

Manual de utilização

Fluent®



Título:	Manual de utiliza	ção do Fluent Dx	Número da peça:	30257931.00	
ID:	403096, pt, V1.0			Traduzido de:	403096, en, V1.0
Versão:	Revisão: Data de Histórico do emissão:			cumento:	
1	0 2025-02-13 Primeira edição				

© 2025, Tecan Trading AG, Switzerland, all rights reserved.

Information contained in this document is subject to change without notice.

TECAN.

Índice

1	Infor	nações sobre este manual	. 7
	1.1	Âmbito deste manual	. 7
	1.2	Operadores económicos	. 7
	1.3	Utilização prevista	. 8
	1.4	Área de utilização/área de aplicação	. 8
	1.5	Utilização inadequada	. 8
	1.6	Garantia	. 9
	1.7	Marcas registadas	. 9
	1.8	Documentos de referência	. 9
	1.9	Cumprimento das leis e das normas	10
	1.10	Convenções do documento	10
-	-		
2	Segu	rança	11
	2.1	Convenções das mensagens de segurança	11
	2.2	Informações gerais de segurança	13
	2.3	Declaração de privacidade da câmara	15
	2.4	Riscos da aplicação	15
	2.5	Empresa exploradora	23
	2.6	Validação do método e do processo	24
	2.7	Qualificações do utilizador	25
	2.8	Elementos de segurança	26
	2.9	Sinais de segurança do produto	32
	2.10	Radiação laser	37
	2.11	Declaração de descontaminação	38
	2.12	Comunicar incidente	39
2	Dede		40
3			4U
	3.1	Placa de características	40
	3.2	Placa do número de série	41
	3.3	Dimensões e pesos	42



	3.4	Fonte de alimentação	43
	3.5	Dados e ligações elétricas	44
	3.6	Condições ambientais	45
	3.7	Emissão e imunidade	46
	3.8	Critérios de aceitação da precisão e exatidão da pipetagem	47
4	Desc	rição do funcionamento	63
	4.1	Vista geral	63
	4.2	Mesa de trabalho	63
	4.3	Braços robóticos	67
	4.4	Sistema de líquido (Liquid FCA)	72
	4.5	Opções e dispositivos	73
5	Elem	entos de controlo	. 85
	5.1	Elementos de comando	85
	5.2	Interface do utilizador	86
	5.3	Sinais de erro e estado do instrumento	91
	5.4	LED de estado do Fluent ID	93
6	Func	ionamento	. 94
	6.1	Instruções de segurança deste capítulo	94
	6.2	Modos de funcionamento	95
	6.3	Colocação em funcionamento	96
	6.4	Antes de iniciar um método	103
	6.5	Executar um método	108
	6.6	Operação DeckCheck	120
	6.7	Method Recovery	122
	6.8	Desligar o instrumento	124
7	Syst	em Care	126
	7.1	Descontaminação	126
	7.2	Agentes de limpeza	127
	7.3	Modo de Cuidados do sistema	128



	7.4	Tabelas de cuidados do sistema	130
	7.5	Atividades de Cuidados do sistema	139
8	Reso	lução de problemas	166
	8.1	Instruções de segurança deste capítulo	166
	8.2	Tabelas da resolução de problemas	166
	8.3	Atividades de resolução de problemas	180
9	Emba	alar, desempacotar, transportar, armazenar e eliminar	210
9	Emba 9.1	alar, desempacotar, transportar, armazenar e eliminar Rótulos das embalagens	210 210
9	Emba 9.1 9.2	alar, desempacotar, transportar, armazenar e eliminar Rótulos das embalagens Eliminação	210 210 211
9	Emba 9.1 9.2	alar, desempacotar, transportar, armazenar e eliminar Rótulos das embalagens Eliminação	210 210 211
9 10	Emba 9.1 9.2 Apoid	alar, desempacotar, transportar, armazenar e eliminar Rótulos das embalagens Eliminação	210210211213
9 10	Emba 9.1 9.2 Apoid 10.1	alar, desempacotar, transportar, armazenar e eliminar Rótulos das embalagens Eliminação D ao cliente Contactos	 210 211 213
9 10	Emba 9.1 9.2 Apoid 10.1	alar, desempacotar, transportar, armazenar e eliminar Rótulos das embalagens Eliminação D ao cliente Contactos	 210 211 213

Índice





1 Informações sobre este manual

Este Operating Manual faculta uma descrição abrangente do sistema Dx do Fluent (simplesmente designado como "Fluent" em todo o manual) e inclui todas as informações necessárias para a sua operação segura e manutenção adequada. Este manual deve ser lido atentamente antes de realizar qualquer trabalho no Fluent e antes de o utilizar.

Este capítulo descreve a finalidade deste manual e especifica o produto a que se refere. Além disso, explica a utilização dos símbolos e convenções, bem como outras informações gerais.

Este manual diz respeito ao próprio Fluent. Para obter informações importantes sobre os submódulos, consulte os manuais específicos dos módulos.



Este Operating Manual não contém qualquer descrição de software. Para obter mais informações sobre o software, consulte o manual do software correspondente. Consulte a secção "Documentos de referência" [> 9].

1.1 Âmbito deste manual

Este manual aplica-se a:

- Fluent Dx 480 (número de peça 30042094)
- Fluent Dx 780 (número de peça 30042095)
- Fluent Dx 1080 (número de peça 30042096)

1.2 Operadores económicos

1.2.1 Fabricante

Endereço do fabricante



Tecan Schweiz AG Seestrasse 103 CH-8708 Männedorf Suíça

1.2.2 Representante Europeu Autorizado

Endereço do Representante Europeu Autorizado



Tecan Austria GmbH

Untersbergstrasse 1a A-5082 Grödig Áustria

1.2.3 Representante Autorizado no Reino Unido

Endereço do Representante Autorizado no Reino Unido

Tecan UK Ltd.

Theale Court 11-13 High Street Theale, Reading, RG7 5AH Reino Unido



1.2.4 Promotor TGA australiano

Endereço do promotor TGA australiano

Promotor australiano Emergo Australia Level 20, Tower II Darling Park 201 Sussex Street Sydney, NSW 2000 Austrália

1.3 Utilização prevista

O Fluent é uma plataforma laboratorial de manuseio de líquidos automatizada para utilização em diagnóstico in vitro. O produto destina-se à automatização da preparação de amostras clínicas e do processamento de ensaios de diagnóstico clínico utilizando amostras humanas. O tipo de amostra e o protocolo de diagnóstico específico são definidos e validados pelo utilizador para o ensaio clínico selecionado. O produto destina-se a ser utilizado por profissionais de saúde e em laboratórios profissionais por pessoal com formação. O produto não se destina a autotestes nem a testes realizados junto do paciente.

Informações adicionais relativas à utilização prevista

Se uma opção ou dispositivo para utilização apenas em investigação (RUO) estiver integrado com o Fluent, a utilização prevista muda para apenas para utilização em investigação. Não deve ser utilizado em procedimentos de diagnóstico.

Se o MCA 384 estiver integrado, a utilização prevista muda para utilização geral (GP).

1.4 Área de utilização/área de aplicação

O Fluent pode ser utilizado em diversos ambientes laboratoriais, de acordo com a sua utilização prevista.

Em cada ambiente, o próprio laboratório é responsável pela validação do instrumento Fluent em conjunto com os líquidos e materiais laboratoriais específicos utilizados no método ou curso de trabalho da aplicação do laboratório.

1.5 Utilização inadequada

A utilização inadequada pode prejudicar o conceito de segurança do Fluent.

- O Fluent não deve ser utilizado com opções ou componentes não aprovados pela Tecan.
- O Fluent não é à prova de explosão e não deve ser instalado em locais onde exista o perigo de explosão.



• O Fluent não deve ser utilizado na ausência de dispositivos de segurança funcionais.

1.6 Garantia

O Fluent não deve ser utilizado com opções ou componentes não aprovados pela Tecan.

A utilização de componentes não aprovados pode comprometer o conceito de segurança do Fluent.

A utilização de componentes não aprovados invalida qualquer garantia de segurança e o cumprimento das normas nacionais e internacionais, conforme exigido pela certificação NRTL, bem como pelas diretivas da CE, entre outros.

1.7 Marcas registadas

Os nomes de produtos, quer sejam marcas registadas ou não registadas, mencionados neste manual são reproduzidos apenas para fins de identificação e permanecem na propriedade exclusiva dos seus respetivos proprietários. Por motivos de simplicidade, os símbolos de marca registada como [®] e [™] não são repetidos no manual.

1.8 Documentos de referência

Esta secção faculta uma lista dos documentos que são necessários ou poderão ser úteis durante a utilização do Fluent.

As ID de documentos listadas abaixo são números de raiz. Consequentemente, não contêm informação sobre a língua, versão do documento ou forma do documento (forma de armazenamento de dados, impressão, ficheiro de download, etc.).



Na base da configuração da sua encomenda, os manuais de utilização para os equipamentos opcionais também se aplicam.

Verifique o âmbito do respetivo documento para se certificar de que possui a versão correta.

A ID do documento não se refere a informações de encomenda. Ao realizar encomendas, consulte o número na capa, caixa do CD, etc.

1.8.1 Manuais do instrumento

- Manual de utilização do Fluent[®] Dx (ID doc. 403096)
- Manual de referência do Fluent[®] Dx (ID doc. 403190)

1.8.2 Manuais de software

- Manual do software adicional de monitorização de amostras Tecan (ID doc. 393933)
- Manual do software de aplicação FluentControl (ID doc. 399935)
- Manual do software de aplicação Introspect (ID doc. 400733)
- Manual do software de aplicação MissionControl (ID doc. 401940)
- Manual de utilização segura do Fluent (ID doc. 403097)



1.8.3 Manuais dos kits de CQ

- Manual de aplicação do kit de CQ (ID Doc. 397069)
- Manual de software de aplicação do kit de CQ (ID Doc. 397070)

1.8.4 Outros documentos de referência

- Manual de utilização do tapete rolante Fluent[®] (ID Doc. 398350)
- Filtro HEPA (ID Doc. Caron 70072)
- Manual de aplicação do Frida Reader™ (ID Doc. 401882)
- Manual de utilização do Te-Shake™ (ID Doc. 391496)
- Manual de utilização do Te-VacS™ (ID Doc. 391236)
- Manual de utilização do Fluent[®] Stacker (ID Doc. 398658)
- Manual de utilização do MIO2 (ID Doc. 394934)
- Manual de utilização do Resolvex i300 (ID Doc. 402756)

1.9 Cumprimento das leis e das normas

As declarações e certificações seguintes aplicam-se ao Fluent:

- Declaração de conformidade CE com as diretivas da UE aplicáveis (marcação CE)
- Certificação de laboratório de testes reconhecido a nível nacional (NRTL)
- Certificação ao abrigo do programa (IECEE) CB (marca CB)

Para obter mais informações sobre a marcação, consultar a secção Placa de características.

1.10 Convenções do documento

Referências As referências cruzadas são apresentadas do seguinte modo - por ex.: cruzadas Consulte a secção "Segurança" [▶ 11] "Segurança" refere-se ao cabeçalho da secção correspondente O número da página é indicado entre parêntesis retos • Pré-requisitos Os pré-requisitos são apresentados do seguinte modo - por ex.: ✓ As "Informações gerais de segurança" foram lidas. **Sugestões** As sugestões adicionais são apresentadas do seguinte modo - por ex.: Para mais informações sobre as convenções e símbolos de segurança, consulte o capítulo "Segurança" [> 11]. llustrações As ilustrações podem apresentar versões dos componentes que não são relevantes para o seu Fluent.



2 Segurança

Este capítulo descreve o conceito de segurança do Fluent, faculta as regras gerais de comportamento correto e adverte sobre os riscos associados à utilização do Fluent.

2.1 Convenções das mensagens de segurança

2.1.1 Palavras-sinal

Tab. 1: Palavras-sinal

Palavra-sinal	Significado
A PERIGO	Indica uma situação perigosa que, se não for evitada, irá causar a morte ou lesões graves.
▲ ATENÇÃO	Indica uma situação perigosa que, se não for evitada, poderá causar a morte ou lesões graves.
▲ CUIDADO	Indica uma situação perigosa que, se não for evitada, poderá causar lesões ligeiras ou moderadas.
AVISO	Indica uma situação que não envolve perigos mas que, se não for evitada, poderá causar danos ou mau funcionamento do equipamento ou resultados de processamento incorretos.



2.1.2 Símbolos de segurança





Campo magnético



2.2 Informações gerais de segurança

ATENÇÃO

O Fluent foi concebido e construído de acordo com o estado da técnica atual e os regulamentos técnicos de segurança reconhecidos. No entanto, podem surgir riscos para os utilizadores, para a propriedade e para o ambiente caso o Fluent seja utilizado sem o devido cuidado e atenção.

A segurança de todos os utilizadores e do pessoal depende do cumprimento rigoroso das instruções de segurança e da consciencialização sobre as advertências relacionadas com a segurança disponibilizadas neste manual.

- Preste especial atenção às seguintes informações gerais de segurança.
- Este manual deve estar sempre disponível para todas as pessoas que desempenham as tarefas nele descritas.
- Utilize sempre o cabo de alimentação fornecido com o instrumento.
- Não utilize o cabo de alimentação com outros produtos.
- Os regulamentos legais, como as leis locais, estatais e federais relativas à utilização ou aplicação, bem como ao manuseamento, de materiais perigosos relacionados com o Fluent têm de ser rigorosamente cumpridos.
- A empresa exploradora é responsável por definir instruções de acordo com os procedimentos da empresa e com os requisitos legais locais. As instruções disponibilizadas pela empresa exploradora têm de ser rigorosamente cumpridas.
- Respeite as condições ambientais corretas para o armazenamento e funcionamento.
- É proibido efetuar alterações estruturais aos dispositivos de segurança.
- Os dispositivos de segurança danificados devem ser substituídos imediatamente, conforme descrito neste manual.
- O Fluent não pode ser modificado de forma alguma sem a consulta prévia e aprovação por escrito da Tecan. As modificações autorizadas ao sistema apenas poderão ser realizadas por um FSE certificado para a reparação e atualização do Fluent. A Tecan rejeitará qualquer reclamação decorrente de modificações não autorizadas.
- Perigo de incêndio causado pela utilização inadequada do Fluent. O Fluent não deve ser instalado em locais onde exista o perigo de explosão.
- Perigo de incêndio causado por líquidos inflamáveis ou pelo líquido do sistema.
- Evite a formação e a acumulação de vapores inflamáveis.
- Às substâncias usadas ou às amostras e reagentes processados no Fluent podem estar associados perigos químicos, biológicos e radioativos (por exemplo, ao carregar e descarregar). O mesmo se aplica à eliminação de resíduos.
 - Tenha sempre atenção aos possíveis perigos associados a estas substâncias.
 - Use vestuário, óculos, respiradores e luvas de proteção adequados.



- A manipulação de substâncias e a eliminação de resíduos poderão estar sujeitos às leis locais, estatais ou federais ou aos regulamentos relativos à saúde, ambiente ou segurança. Cumpra rigorosamente as disposições correspondentes.
- Todas as contaminações têm de ser tratadas imediatamente conforme descrito neste manual.
- O utilizador é responsável por assegurar que o Fluent é sempre utilizado nas condições adequadas e que a manutenção, a assistência e as reparações são sempre realizadas com cuidado, atempadamente e apenas por pessoal autorizado.
- Risco de resultados de medição incorretos. Depois da realização dos cuidados do sistema ou da manutenção, o funcionamento só deve ser retomado depois de terem sido verificadas as corretas condições de utilização do sistema.
- Para assegurar um bom desempenho do sistema e a sua fiabilidade, na manutenção e reparação, utilize sempre consumíveis recomendados dentro do prazo de validade e peças sobresselentes originais.
- Se a pele entrar em contacto com o líquido do sistema no instrumento, poderão ocorrer lesões.
 - Usar sempre vestuário de proteção, de acordo com as BPL.
- Carga pesada! Não levantar o instrumento.
- Não operar o sistema sem os tabuleiros de recolha e os segmentos da plataforma.
- Os tabuleiros de recolha recolhem derrames de líquidos que podem ocorrer na área de carregamento manual da plataforma. O sistema deve ser operado com o maior número possível de tabuleiros de recolha instalados sob a plataforma para recolher todos os derrames de líquidos. Não operar o sistema sem os tabuleiros de recolha.
- Se não forem tolerados resíduos de transição, recomendamos vivamente a utilização de pontas descartáveis com filtro.
- Possível colisão. Não colocar dispositivos na plataforma sem dados do modelo Tecan.
- A extensão 300 destina-se a uma carga máxima de 40 kg (88 lbs) e apenas à utilização com opções com menos de 40 kg (88 lbs).
- O Fluent é fornecido com um sinal de segurança de perigo biológico que deve ser aplicado pelo utilizador em caso de utilização de substâncias biológicas perigosas. Aplicar a etiqueta na porta dianteira numa posição visível para o utilizador e conveniente para a aplicação. Consulte a secção "Sinais de segurança do produto" [▶ 32].
- As opções utilizadas na mesa de trabalho do Fluent podem gerar campos magnéticos fortes, que podem interferir com o funcionamento de dispositivos médicos implantados ou utilizados por um operador, como pacemakers ou bombas de insulina. O Fluent é fornecido com um sinal de segurança de campos magnéticos fortes que deve ser aplicado pelo utilizador na porta frontal, numa posição visível para o utilizador e adequada para a aplicação em caso de utilização de opções que gerem campos magnéticos fortes.
- O cabo Ethernet das câmaras DeckCheck será instalado por um FSE e tem de estar sempre instalado no PC com FluentControl (CEM). A interface Ethernet não pode ser ligada a uma rede.



2.3 Declaração de privacidade da câmara

O sistema Fluent está equipado com câmaras montadas no interior da parte da frente do equipamento. As câmaras estão focadas na mesa de trabalho de na mesa de trabalho posterior. É possível obter vistas para baixo através dos painéis laterais de vidro acrílico.

- O utilizador é responsável por informar as pessoas presentes na divisão de que as câmaras estão em funcionamento.
- O utilizador é responsável por garantir que não é possível identificar os técnicos nas imagens capturadas, por exemplo, se o instrumento estiver adjacente (de lado) a uma secretária ou se forem efetuados recortes nos painéis laterais ou traseiro, ou ainda se for utilizado um painel de vidro acrílico para substituir a parede traseira.

Função do sistema/ módulo	Modo de falha possível	Potencial efeito da falha	Causa possível/ potencial	Identificação ou atenuação
Sistema	Manutenção insuficiente	Segurança ou saúde dos utilizadores: possível contaminação do instrumento	Erro de utilização: ignorar o manual de utilização ou as instruções de manutenção	O utilizador tem de garantir a utilização de consumíveis adequados e o cumprimento das instruções de manutenção preventiva (consulte a secção "System Care" [▶ 126]). O utilizador deve utilizar vestuário,
				luvas e óculos de proteção de acordo com as BPL e os regulamentos locais em vigor.
Sistema	Incêndio	Segurança ou saúde dos utilizadores: incêndio no laboratório do operador (incineração de instrumento)	Gás dos líquidos inflamáveis voláteis; faíscas irradiadas pelo quadro elétrico	O instrumento não é à prova de explosão e o cliente deve certificar-se de que não existe uma concentração elevada de vapor (consulte a secção "Informações gerais de segurança" [▶ 13]).
Módulo FCA e Air FCA	Desgaste dos mecanismos do eixo Z (utilização acima da média)	Segurança ou condições clínicas da amostra: posicionamento Z potencialmente errado no material laboratorial	Utilização do dispositivo acima da média em combinação com a utilização de pontas descartáveis Elevada percentagem de passos de perfuração na aplicação	O sistema informa o utilizador se os eixos Z tiverem atingido 90% da vida útil esperada.

2.4 Riscos da aplicação



Função do sistema/ módulo	Modo de falha possível	Potencial efeito da falha	Causa possível/ potencial	Identificação ou atenuação
Módulo FCA e Air FCA	Desgaste dos mecanismos do eixo P (utilização acima da média)	Segurança ou condições clínicas da amostra: posicionamento P potencialmente errado no material laboratorial	Utilização do dispositivo acima da média em combinação com a utilização de pontas descartáveis Elevada percentagem de passos de perfuração na aplicação	O sistema informa o utilizador se os eixos P tiverem atingido 90% da vida útil esperada.
Módulo FCA e Air FCA	Abrasão da roda dentada do acionamento X (utilização acima da média)	Segurança ou condições clínicas da amostra: potencial contaminação das amostras com partículas de poliamida	Utilização do dispositivo acima da média em combinação com a colocação de material laboratorial crítico na parte posterior do instrumento	Evitar colocar elementos sensíveis a partículas (por exemplo, amostras e reagentes) na traseira do instrumento ou colocar a proteção contra partículas (ou seja, tampas) sobre o material laboratorial.
Módulo FCA e Air FCA	Sinais com interferência devido a perfuração do septo	Segurança ou condições clínicas da amostra do paciente: cLLD errada, dando origem a aspiração de ar e a potenciais resultados falsos	Interação da ponta com o septo/ película	Trabalhar apenas com películas não condutoras em aplicações de perfuração, em conjunto com a deteção do nível de líquido no FCA e no Air FCA. Consultar o manual de referência. O utilizador tem de validar a deteção do líquido em combinação com a perfuração para o FCA e o Air FCA.
Módulo FCA e Air FCA	Tratamento errado da amostra, cLLD errada devido a espuma ou bolhas no frasco de reagente	Segurança do processo: amostras processadas de forma errada	As bolhas ou espuma no frasco de reagente causam uma cLLD errada e uma potencial aspiração de ar com o FCA e o Air FCA	O utilizador é responsável pela validação da aplicação/do processo relativamente à preparação adequada da amostra.



Função do sistema/ módulo	Modo de falha possível	Potencial efeito da falha	Causa possível/ potencial	Identificação ou atenuação
Módulo FCA e Air FCA	Bloqueio da ponta	Segurança ou condições clínicas da amostra do paciente: volume pipetado potencialmente errado	Aspiração no fundo do poço (bloqueio da ponta)	O utilizador tem de validar a aplicação para prevenir a aspiração demasiado próxima do nível máx. Z do material laboratorial personalizado.
Módulo FCA e Air FCA	Sistema de tubagem do FCA: crescimento de microrganis mos	Segurança ou condições clínicas da amostra do paciente: volume pipetado errado ou contaminação das amostras	Crescimento de microrganismos (biofilme na superfície interior)	Utilize água desionizada como líquido do sistema para o FCA e execute a manutenção diária para lavar o sistema de acordo com as instruções de cuidados diários do sistema (consulte a secção "System Care" [▶ 126]), respeitando também os agentes de limpeza permitidos para esse passo.
Módulo MCA 96	Extravasame nto do líquido da amostra na microplaca durante a pipetagem	Segurança ou condição clínica da amostra do paciente: potencial contaminação cruzada das amostras (extravasament o)	Níveis Z definidos de forma errada pelo utilizador (por exemplo, aspiração a partir da posição máx. Z)	Definir posições seguras para aspiração e distribuição. Consultar o manual de referência.
Módulo MCA 96	Amostras falham completa ou parcialmente a posição prevista no modo de distribuição livre	Segurança ou condição clínica da amostra do paciente: potencial contaminação cruzada	Cargas eletrostáticas na extremidade da ponta devido à utilização do instrumento fora das condições especificadas levam a que a amostra permaneça suspensa na ponta ou a pulverizações não controladas	O utilizador tem de respeitar as condições de funcionamento especificadas para o manuseamento de líquidos do MCA, especialmente as instruções sobre a humidade mínima necessária (consulte a secção "Condições ambientais" [▶ 45]). O utilizador tem de definir a altura de distribuição sempre dentro do poço. Consultar o manual de referência.



Função do sistema/ módulo	Modo de falha possível	Potencial efeito da falha	Causa possível/ potencial	Identificação ou atenuação
Módulo MCA 96	Misturar ar em vez de líquido (amostra/ reagente) para a pipetagem da mistura	Segurança ou condição clínica da amostra do paciente: amostras processadas de forma potencialmente incorreta, causando resultados falsos	Parâmetros de monitorização inadequados devido à combinação errada de pontas e microplacas	O utilizador deve comparar a mesa de trabalho real e virtual, ao utilizar o nome do material laboratorial na mesa de trabalho virtual. O utilizador deve observar o design de cor única (específica para o tipo de ponta) e a identificação (para filtro e não filtro) das caixas DiTi. O utilizador tem de verificar a disposição da mesa de trabalho antes de iniciar um processo.
Específico para ponta descartável	Obter DiTi: está montado o tipo de ponta incorreto	Segurança ou condição clínica da amostra do paciente: Possibilidade de aspiração ou aspiração insuficiente da amostra Potencial contaminação cruzada das amostras	Erro de utilização: Disposição incorreta da plataforma: o utilizador coloca a caixa de pontas na posição errada: as pontas são menores do que o esperado Disposição incorreta da plataforma: o utilizador coloca na plataforma a caixa de pontas com pontas não filtradas em vez de pontas filtradas Disposição incorreta da plataforma: o utilizador coloca a caixa de pontas na posição errada: a ponta tem um volume inferior ao esperado (por exemplo, 100 µl em vez de 200 µl); comprimento da ponta conforme esperado; líquido aspirado na MCH	O utilizador deve comparar a mesa de trabalho real e virtual, ao utilizar o nome do material laboratorial na mesa de trabalho virtual. O utilizador deve observar o design de cor única (específica para o tipo de ponta) e a identificação (para filtro e não filtro) das caixas DiTi. O utilizador tem de verificar a disposição da mesa de trabalho antes de iniciar um processo. O design mecânico assegura a visibilidade do filtro branco. O manual de referência contém informações sobre o código de cores das caixas DiTi, a diferença de comprimento e as DiTi filtradas. Consultar o manual de referência.



Função do sistema/ módulo	Modo de falha possível	Potencial efeito da falha	Causa possível/ potencial	Identificação ou atenuação
Específico para ponta descartável	Descida incompleta das pontas: algumas pontas contaminada s permanecem suspensas na cabeça e caem nas placas de amostras	Segurança ou condição clínica da amostra do paciente: potencial contaminação cruzada	Cargas eletrostáticas causadas	O utilizador tem de respeitar as condições de funcionamento especificadas para o manuseamento de líquidos do MCA, especialmente as instruções sobre a humidade mínima necessária (consulte a secção "Condições ambientais" [▶ 45]). As pontas descartáveis não se destinam à reutilização.
Módulo RGA	Perda da placa devido a colisão com material laboratorial desalinhado	Segurança do processo: perda da placa, perda das amostras	Se estiverem empilhadas mais de 4 microplacas, pode ocorrer um desalinhamento durante o transporte	Os movimentos da placa têm de ser validados antes de executar scripts com amostras reais. Consultar a lista de verificação de validação no manual do software de aplicação.
Módulo de software FluentContro I	Base da mesa de trabalho: estado da DiTi errado comunicado	Segurança do processo: contaminação cruzada/ resultados errados	Contaminação cruzada devido a informações erradas sobre o estado de utilização das pontas	Não utilizar "Set Tips Back" (Repor pontas) se o modo de falha resultar num risco de gravidade elevada.
Módulo de software FluentContro I	Core.Scriptin g.Programmi ng SetVariable no tempo de execução: valor errado	Segurança do processo: resultados errados	Erro no software: a variável está configurada para um valor errado	Validar a aplicação para a fonte, o destino e os âmbitos da variável específica. Consultar a lista de verificação de validação no manual do software de aplicação.
Módulo de software FluentContro I	Core.Scriptin g.Programmi ng QueryVariabl e no tempo de execução ou no início do script: erro de apresentaçã o da IU/ aceitação do valor da IU	Segurança do processo: resultados errados	O valor numérico está formatado ou convertido de forma errada na IU	Validar a aplicação para a fonte, o destino e os âmbitos da variável específica. Consultar a lista de verificação de validação no manual do software de aplicação.



Função do sistema/ módulo	Modo de falha possível	Potencial efeito da falha	Causa possível/ potencial	Identificação ou atenuação
Módulo de software FluentContro I	Core.Scriptin g.Programmi ng ImportVariabl e no tempo de execução: valor errado importado	Segurança do processo: resultados errados	O valor errado é recuperado da fonte da importação	Validar a aplicação para a fonte, o destino e os âmbitos da variável específica. Consultar a lista de verificação de validação no manual do software de aplicação.
Módulo de software FluentContro I	Core.Scriptin g.Programmi ng ExportVariab le no tempo de execução: valor errado exportado para o ficheiro	Segurança do processo: resultados errados	O valor errado está gravado no ficheiro de exportação	Validar a aplicação para a fonte, o destino e os âmbitos da variável específica. Consultar a lista de verificação de validação no manual do software de aplicação.
Módulo de software FluentContro I	API: obter/ definir a variável ou resolver as falhas da expressão	Segurança do processo: resultados errados	Valor da variável errado recuperado ou valor errado atribuído/resultado da expressão errado enviado	Validar a aplicação para a fonte, o destino e os âmbitos da variável específica. Consultar a lista de verificação de validação no manual do software de aplicação.
Luz UVC	Utilização errada na aplicação	Falta de eficácia	Utilização errada na aplicação	Consulte as instruções específicas na secção "Radiação ótica (UVC)" [▶ 31].
Rotador de tubos/Pontas de perfuração (mistura e perfuração)	Utilização errada na aplicação	Falta de eficácia	Utilização errada na aplicação	Consulte as instruções específicas na secção "Mistura e perfuração" [▶ 77].
Leitor Frida	Utilização errada na aplicação	Falta de eficácia	Utilização errada na aplicação	Consulte as instruções específicas na secção "Frida Reader" [▶ 80].
Qualquer	Utilização ineficaz na aplicação	Falta de eficácia na aplicação	Falta de cuidados do sistema	Consulte as instruções específicas do componente no capítulo "System Care" [▶ 126]
Processame nto de materiais potencialmen te perigosos	Contaminaçã o com materiais potencialmen te perigosos	Riscos potenciais para utilizadores, propriedade e ambiente.	Falta de adesão às informações gerais de segurança	Consulte as instruções específicas do componente na secção "Informações gerais de segurança" [▶ 13].



Função do sistema/ módulo	Modo de falha possível	Potencial efeito da falha	Causa possível/ potencial	Identificação ou atenuação
MCA 96, cLLD	Resultado de medição incorreto	Medição errada: o nível de líquido detetado não se aplica a todos os poços da microplaca: resultado de teste errado ou perda de amostra	Utilização de material laboratorial inadequado/ incompatível	A função de cLLD só pode ser utilizada em recipientes.
MCA 96, geral	Incompatibili dade de consumíveis/ componente s/módulos	Perda de DiTi durante o processo: potencial perda irrecuperável de amostras. Potencial contaminação cruzada.	Utilização de consumíveis inadequados/ incompatíveis. É comandada uma recolha de compensação que não é compatível com a caixa ou tabuleiro a partir do qual as DiTi devem ser montadas (por exemplo, versão errada do tabuleiro das DiTi). As DiTi estão montadas incorretamente e perdem-se durante o processo	Utilizar caixas DiTi com tabuleiro combinado Tecan.
MCA 96, cLLD	Resultado de medição incorreto	Medição errada: deteção de falsos positivos do subsistema de cLLD.	O utilizador coloca na mesa de trabalho DiTi diferentes das refletidas no software. O número de DiTi montadas difere do número de DiTi esperadas pelo software devido a uma ação do utilizador (por exemplo, o utilizador remove algumas DiTi). Deteções de falsos positivos devido a um limiar de cLLD incorreto.	Importante: o número de DiTi utilizadas para a deteção de cLLD tem de corresponder ao número refletido no software.



Função do sistema/ módulo	Modo de falha possível	Potencial efeito da falha	Causa possível/ potencial	Identificação ou atenuação
MCA 96, geral	Contaminaçã o da amostra	Contaminação da amostra devido a derrame após uma colisão. Resultado de teste errado ou perda de amostra.	Parâmetros errados tomados (vetor errado para a recolha ou entrega, tipo de ferramenta errado). Colisão da cabeça do MCA 96 durante o movimento do vetor. Por exemplo, com material laboratorial elevado em estações adjacentes.	Existem riscos associados à utilização incorreta da funcionalidade de movimentação do vetor.
MCA 96, geral	Contaminaçã o da amostra	Segurança ou condição clínica da amostra do paciente: Contaminação da amostra: potencial contaminação cruzada devido a abrasão da correia do manipulador.	Abrasão do material da correia ao longo da vida útil: as partículas de abrasão da correia do eixo G podem cair para as amostras de pacientes na mesa de trabalho e acabar por contaminar a amostra quimicamente.	Partículas de abrasão (das correias) e pó podem cair para a mesa de trabalho e contaminar as amostras/os produtos químicos. Para evitar esta situação, o material laboratorial que contém tais líquidos sensíveis deve ser protegido através da colocação de tampas.
MCA 96, geral	Contaminaçã o da amostra	Segurança ou condição clínica da amostra do paciente: Resultados potencialmente errados: potencial contaminação cruzada resultante de derrames para cavidades adjacentes da placa.	Salpicos/Derrame de líquido da amostra. Contaminação dos cones de DiTi por líquido no material laboratorial no manipulador, por exemplo, quando um material laboratorial cheio é colocado no lixo e ocorre o derrame de líquidos.	Esvaziar o material laboratorial antes de o eliminar.



Função do sistema/ módulo	Modo de falha possível	Potencial efeito da falha	Causa possível/ potencial	Identificação ou atenuação
MCA 96, geral	Transferênci a de amostras ou reagentes (FC)	Cones de DiTi e bloco de cilindros contaminados, resultando na potencial contaminação cruzada das amostras.	O dispositivo tenta recolher as DiTi. Uma vez que o software não consegue detetar DiTi em falta, este continua sem erros. Os cones de DiTi podem entrar em contacto e aspirar o líquido do recipiente	O MCA 96 não consegue detetar caixas de pontas vazias ao utilizar tipos de caixas de pontas em que a extremidade da ponta esteja nivelada com a superfície superior da caixa.
MCA 96, geral	Salpicos/ Derrame de líquido da amostra	Contaminação da amostra devido a derrame por falta de material laboratorial suficientemente rígido.	Salpicos/Derrame de líquido da amostra devido à utilização de material laboratorial não suficientemente rígido.	Para evitar qualquer deformação do material laboratorial, utilizar apenas material laboratorial suficientemente rígido para transportar líquidos.
MCA 96 geral	O dedo do manipulador pode ficar danificado após uma colisão	Dedo do manipulador danificado		Verificar os dedos do manipulador e, se estes apresentarem danos, substituir.
Mistura e perfuração da estação de lavagem	Contaminaçã o da amostra	Os resíduos não estavam a ser drenados corretamente. Contaminação da amostra devido a drenagem incorreta dos resíduos e lavagem incorreta das pontas.	Pontas incorretamente lavadas devido a extravasamento da estação de lavagem causado por obstrução da ligação à estação de lavagem.	O sistema deve ser submetido a manutenção regular. Os conectores da estação de lavagem devem ser trocados em ciclos de 2-3 meses para aplicações de sangue total de modo a evitar obstruções. A drenagem de líquidos potencialmente corrosivos, como lixívia a 2%, através da estação de lavagem, dos conectores e dos tubos não deve ser efetuada sem lavagem adicional com líquidos neutros, como água, para evitar a corrosão

2.5 Empresa exploradora

A empresa exploradora deve garantir que o Fluent e, em especial, as características de segurança, funcionam adequadamente e que todo o pessoal em contacto com o instrumento recebe formação adequada.



Responsabilida des

- Validação do método e do processo.
- Definir os processos em conformidade com procedimentos operativos normalizados.
- Garantir que as qualificações de instalação e operacionais (QI QO) foram realizadas.
- Garantir que todo o pessoal em contacto com o Fluent recebe formação adequada.
- Garantir que está disponível vestuário e equipamento de proteção adequado.
- · Garantir a manutenção e o funcionamento seguro do Fluent.
- Exigir o cumprimento dos regulamentos e diretivas de segurança laboratorial.

2.6 Validação do método e do processo

Ao realizar a validação do método e do processo, ter em atenção o seguinte:

- Se utilizar pontas fixas com o FCA, certifique-se de que o procedimento de lavagem é eficaz para o intervalo de concentração esperado da amostra e para a sensibilidade da análise.
- Verificar se os volumes pipetados cumprem os requisitos de precisão e exatidão do processo a ser automatizado.
- Ao utilizar material laboratorial personalizado ou que não seja da Tecan e ao aspirar com monitorização, garantir que a definição do recipiente é a correta (ou seja, que é utilizada a velocidade correta para a monitorização) para evitar a aspiração de ar.
- A funcionalidade do separador de fases foi verificada para utilização com pontas descartáveis Tecan padrão de 1 ml e pontas descartáveis Tecan de abertura larga de 1 ml. Para obter mais informações sobre os consumíveis Tecan suportados, consulte o manual de referência (consulte a secção "Documentos de referência" [> 9]).
- Validar a deteção de líquido na estação de transferência Fluent Stacker.
- Validar a utilização correta da estação de lavagem do MCA pela aplicação.
- Validar a aplicação em matéria de volumes de pipetagem e monitorização corretos.
- Validar a aplicação para impedir a aspiração demasiado próxima do máx. Z do material laboratorial personalizado.
- Validar as aplicações de perfuração relativamente aos fixadores necessários (ativos ou passivos).
- Se os produtos químicos e o material laboratorial não forem removidos, será necessário avaliar o efeito da luz UVC nos respetivos produtos e materiais presentes na plataforma e validar a análise.
- Incluir uma verificação manual pós-execução para obter os volumes de pipetagem corretos.
- Os técnicos devem estar informados acerca da declaração de privacidade da câmara (consulte a secção "Declaração de privacidade da câmara" [▶ 15]).

Responsabilida de do operador principal



2.7 Qualificações do utilizador

O pessoal de laboratório tem de possuir as qualificações e formação adequadas para operar o Fluent. O trabalho descrito neste Operating Manual só deve ser realizado por pessoal autorizado com as qualificações definidas abaixo.

O pessoal de laboratório deve:

- ter formação técnica adequada,
- · estar familiarizado com os regulamentos e diretivas de segurança laboratorial,
- estar familiarizado com as instruções dos elementos de segurança do instrumento,
- usar vestuário e equipamento de proteção,
- estar familiarizado com as boas práticas laboratoriais e cumpri-las
- e ter lido e compreendido as instruções do manual de utilização.

A Tecan recomenda que o operador frequente um curso de formação para operadores. Peça à Assistência a clientes da Tecan informações sobre os cursos disponíveis. Consulte a secção "Apoio ao cliente" [> 213].

2.7.1 Operador

O operador (técnico laboratorial) trabalha para a empresa exploradora.

- Nenhum conhecimento específico sobre a aplicação ou o sistema
- Domínio das línguas locais
- O domínio do inglês é preferencial

O operador tem direitos de acesso ao software da aplicação que lhe permitem executar métodos e efetuar os cuidados do sistema e receberá a formação necessária ministrada pelo operador principal.

2.7.2 Operador principal

O operador principal (especialista na aplicação) dá assistência à empresa exploradora ou trabalha para a mesma empresa.

 Competências necessárias
 • Vasto conhecimento da aplicação

 • Conhecimento limitado do sistema

 • Domínio das línguas locais

 • Domínio do inglês

 • Conhecimento aprofundado do manual do software correspondente

 • Instruir o operador

Responsabilida des

- Escrever, executar e validar métodos
- Ajudar o operador a resolver problemas com o instrumento

Competências necessárias



2.8 Elementos de segurança

Componentes em movimento

Os elementos de proteção e segurança instalados no Fluent não devem ser removidos, desativados ou substituídos durante o funcionamento.

 Se algum dos dispositivos for removido (por exemplo, para trabalhos de manutenção), todos os dispositivos de proteção e segurança devem ser novamente instalados, reativados e verificados antes de retomar as operações.

Os painéis de segurança e os sensores de segurança são partes integrantes do Fluent, enquanto os bloqueios da porta do instrumento e os bloqueios da porta do armário podem ser incluídos apenas em determinadas configurações do sistema.

2.8.1 Painéis de segurança

O Fluent está protegido com painéis de segurança:

O **painel de segurança frontal** pode ser aberto e está equipado com sensores de porta que desencadeiam uma paragem ativa. O painel de segurança frontal pode ser bloqueado com fechos de porta opcionais.

O Fluent com um braço MCA 96, pontas de perfuração ou Resolvex i300 só pode ser utilizado com o painel de segurança frontal completo.

A abertura da porta é apoiada por molas a gás. Para uma segurança ideal e acesso total ao instrumento, o operador tem de abrir completamente a porta antes de continuar.

O **painel do diluidor** pode ser aberto sem afetar o funcionamento do Fluent (exceto para Fluents com a opção de luz UVC instalada – o sensor do painel do diluidor irá acionar uma paragem rápida quando o painel for aberto).

Os painéis de segurança superior e lateral estão fixos.

2.8.1.1 Painéis de segurança frontais

O painel de segurança frontal impede o acesso direto aos braços robóticos e aos elementos da mesa de trabalho do instrumento durante o funcionamento. Esta configuração existe para benefício da segurança individual e melhora a segurança do método. Além disso, o painel de segurança frontal protege o utilizador contra o derrame de amostras ou reagentes. Existem diferentes tipos de painéis frontais de segurança.



Painel de segurança frontal total



Fig. 1: Painel de segurança frontal total

O painel de segurança frontal total possui as seguintes características:

- Sem acesso aos componentes em movimento (componentes em movimento, perigo mecânico)
- Proteção das amostras contra influências exteriores (segurança do método)
- Proteção contra o derrame de amostras ou reagentes



Com os painéis de segurança frontais totais, é possível apenas o carregamento revisto por lote.

2 - Segurança Elementos de segurança





Painel de segurança frontal total (UVC)

Fig. 2: Painel de segurança frontal total (UVC)

O painel de segurança frontal total (UVC) possui as seguintes características:

- Sem acesso aos componentes em movimento (componentes em movimento, perigo mecânico)
- Proteção das amostras contra influências exteriores (segurança do método)
- · Proteção contra o derrame de amostras ou reagentes
- Proteção contra a radiação ótica (UVC)



Com os painéis de segurança frontais totais, é possível apenas o carregamento revisto por lote.

Componentes em movimento!

O MCA, o FCA e o Air FCA em movimento podem causar ferimentos nas mãos ao colocá-las no instrumento, através do meio painel de segurança frontal ou do painel de segurança frontal com expansão, durante uma execução.

Não coloque as mãos dentro do instrumento durante uma execução.



Meio painel de segurança frontal



Fig. 3: Meio painel de segurança frontal

O meio painel de segurança possui as seguintes características:

- Acesso restrito aos componentes em movimento (componentes em movimento, perigo mecânico)
- Proteção contra o derrame de amostras ou reagentes



Com o meio painel de segurança frontal, o operador tem acesso restrito à mesa de trabalho do instrumento. É possível carregar e descarregar as calhas sem abrir o painel, ou seja, o operador tem a possibilidade de voltar a carregar amostras ou reagentes durante a execução do método.

2 - Segurança Elementos de segurança



Painel de segurança frontal com expansão



Fig. 4: Painel de segurança frontal com expansão

O painel de segurança possui as seguintes características:

- Acesso restrito aos componentes em movimento (componentes em movimento, perigo mecânico)
- Proteção contra o derrame de amostras ou reagentes
- Permite a utilização de uma estação de resíduos DiTi frontal, que sobressai da mesa de trabalho e requer uma abertura voltada para baixo no painel de segurança frontal.



Com o painel de segurança frontal com expansão, é possível apenas o carregamento revisto por lote.

2.8.1.2 Painéis de segurança para dispositivos opcionais

Caso um dispositivo opcional seja adicionado ou removido da lateral do Fluent, é necessário instalar um painel de segurança lateral adequado. Consulte o "Apoio ao cliente" [> 213].

2.8.2 Bloqueios da porta do instrumento (opcional)

Dois bloqueios opcionais da porta do instrumento podem impedir a abertura do painel de segurança e protegem o processo em curso. Tal previne a interrupção injustificada da execução do processo. Para parar um processo, pode introduzir um pedido de pausa através do ecrã tátil.



Determinadas configurações requerem fechos das portas. Se uma configuração for alterada, poderá ser necessária uma atualização com fechos das portas. Consulte a secção "Apoio ao cliente" [> 213].

2.8.3 Bloqueios da porta do armário

Se um eixo longo do RGA tiver acesso abaixo da plataforma, a porta do armário mais próxima do ponto de acesso deve estar equipada com a opção de um sensor de bloqueio da porta. Caso estejam implementados mais do que um ponto de acesso por baixo da plataforma ou se o ponto de acesso for alterado ao longo da existência do instrumento, cada porta junto do ponto de acesso deve estar equipada com um sensor de bloqueio da porta.

Se o instrumento tiver uma tampa HEPA ou um Resolvex i300, todas as portas do armário devem estar equipadas com um sensor de fecho da porta.

2.8.4 Radiação ótica (UVC)

O Fluent pode ser equipado com um exaustor com filtro HEPA opcional, que inclui uma luz UVC ou uma opção de luz UVC separada.

A exposição à radiação da luz UVC tem de ser evitada, uma vez que esta pode provocar lesões. A luz UVC desliga-se automaticamente quando o painel de segurança frontal é aberto e, no caso da opção de luz UVC, também quando a tampa do diluidor é aberta. Estão instalados no Fluent painéis de segurança especiais resistentes aos raios UVC, em conjunto com a luz UVC.

A luz UVC pode ser utilizada em procedimentos de descontaminação. A adequação e eficácia da utilização de raios UVC para processos individuais tem de ser validada pelo utilizador.



Consultar também o manual disponibilizado pelo fabricante do exaustor com filtro HEPA.

2.8.5 Bloqueios da porta externos

Os bloqueios da porta externos serão implementados nas instalações do Fluent num invólucro externo. Os painéis da porta do invólucro externo substituem a função de segurança mecânica do painel de segurança frontal e das portas do armário do Fluent e todos os bloqueios da porta externos com sensores integrais substituem o sensor da porta e as funções de bloqueio da porta do painel de segurança frontal e das portas do armário do Fluent.



Os bloqueios da porta externos não permitem uma paragem ativa. Para parar ou colocar em pausa o processo, pode introduzir um pedido de pausa através do ecrã tátil.



2.9 Sinais de segurança do produto

Os sinais de segurança estão afixados no Fluent por motivos de segurança. Sinais de segurança danificados, em falta ou ilegíveis têm de ser substituídos imediatamente, conforme ilustrado. Para conhecer o significado dos símbolos de segurança, consulte a secção "Convenções das mensagens de segurança" [▶ 11].



UVC



Fig. 5: Instrumento padrão



Fig. 6: Instrumento com UVC





Fig. 7: Vista do interior



Fig. 8: Perigo biológico







Instrumento com meio painel de segurança frontal O Fluent é fornecido com um sinal de segurança de perigo biológico que deve ser aplicado pelo utilizador em caso de utilização de substâncias biológicas perigosas.

Aplicar a etiqueta na porta dianteira numa posição visível para o utilizador e conveniente para a aplicação.



Fig. 9: Instrumento com meio painel de segurança frontal





Instrumento com painel de segurança frontal com expansão

Fig. 10: Instrumento com painel de segurança frontal com expansão





Fig. 11: Extensão da plataforma



Placa de

segurança do FCA



2.9.1 Estação de trabalho de mistura e perfuração

Fig. 12: Placa de segurança


2.10 Radiação laser

O Fluent pode ser equipado com leitores de códigos de barras a laser. A radiação laser destes leitores de códigos de barras é um raio de baixa potência, colimado no espectro visível. As classes de laser de cada leitor de códigos de barras e de todo o sistema Fluent são indicadas na etiqueta de segurança de laser afixada no hardware correspondente.

Todos os módulos com laser estão marcados com as etiquetas de segurança laser adequadas.

O instrumento Fluent foi testado e certificado de acordo com as normas CEI 60825-1:2007 e CEI 60825-1:2014.



O Fluent é um produto laser de classe 1, de acordo com a norma IEC 60825-1:2014, que emite radiação laser.

O raio laser pode causar encadeamento, cegueira por ofuscamento e imagens persistentes.

• Não olhe para o raio laser nem para os reflexos deste.

2.10.1 Dispositivos com radiação laser

Pode ser montado no dispositivo um leitor de código de barras autónomo.

Certifique-se de que a etiqueta de segurança está sempre corretamente afixada no leitor de código de barras:

- Etiqueta de radiação laser explicativa (A): Identifica um PRODUTO LASER DE CLASSE 2 de acordo com a norma IEC 60825-1 que possui um leitor de código de barras embutido com um laser de baixa potência visível. Instrui o utilizador para não olhar para o raio nem para o reflexo deste.
- Os lasers de classe 2 são operados apenas quando o sistema está a ser executado e não há qualquer interface para o operador.

Colocação da etiqueta	Explicação
A LSER RADIATION D MOT STARE INTO BEAK Dutput 854W Markength 854W Markeng	Leitor de código de barras autónomo instalado num hotel: etiqueta colocada por baixo do leitor.
LASER RADIATION DO NOT STARE ITO DEM Duise duration 11249 CLASS 2 LASER PRODUCT IEC/EN 60825-1: 2007	Leitor de código de barras autónomo instalado no Robotic Gripper Arm: etiqueta colocada no leitor.



Colocação da etiqueta	Explicação
	Loading ID: etiqueta colocada no lado traseiro da caixa do leitor.
ACID EXPOSES - LASER RADATION S BMITTED FROM THIS APERTURE	Loading ID: etiqueta colocada na lateral da caixa do leitor.

2.11 Declaração de descontaminação

Além dos cuidados regulares do sistema, e de acordo com os regulamentos laboratoriais padrão, o Fluent e os seus componentes e acessórios têm de ser completamente descontaminados nas seguintes circunstâncias:

- Antes da realização de qualquer trabalho de manutenção ou de reparação no Fluent e, em especial, antes da intervenção por parte de um FSE no Fluent
- Em caso de acidentes (por ex., embate, derrame, etc.)
- Antes de devolver o Fluent ou os seus componentes ou acessórios, à Tecan (por ex., para reparação)
- Antes do armazenamento
- Antes da eliminação
- De forma geral, antes de deslocar o Fluent ou os seus componentes da respetiva localização

A descontaminação efetiva de todo o equipamento é da total responsabilidade do proprietário do instrumento.

Antes de qualquer intervenção no Fluent por um FSE e antes de devolver o Fluent ou os seus componentes ou acessórios à Tecan, o proprietário do instrumento tem de preencher e assinar o formulário da Declaração de descontaminação, que confirma que a descontaminação foi realizada de acordo com as orientações das boas práticas de laboratório. Contacte o responsável local pela assistência técnica para obter este formulário e consulte a secção Descontaminação.



A Tecan reserva-se o direito de recusar o processamento de qualquer Fluent ou respetivos componentes ou acessórios que não sejam acompanhados pelo formulário da Declaração de descontaminação.



2.12 Comunicar incidente

Qualquer incidente grave que ocorra relacionado com o dispositivo tem de ser comunicado ao fabricante e à autoridade competente do Estado-Membro onde o utilizador e/ou paciente está estabelecido. Consulte a secção Fabricante para obter o endereço do fabricante.



3 Dados técnicos

3.1 Placa de características



Fig. 13: Placa de características

A placa de características encontra-se na parte traseira do Fluent e contém as seguintes informações:

Dados de identificação	Modelo
	REF: informação de encomenda (número de material e nível de revisão)
	Data de fabrico (AAAAMMDD)
	SN: número de série
Dados técnicos	U, f: tensão de alimentação (volts), frequência (Hertz)
	P: consumo de energia (VA)
	Fusível: especificação do fusível
Dados de endereço	Nome e endereço do fabricante
Dados de	Marca de conformidade
conformidade	UDI: Identificação única do dispositivo O símbolo UDI identifica o suporte de dados na etiqueta.
	EC REP: Representante Europeu Autorizado
	IVD: Dispositivo médicos para diagnóstico in vitro



3.2 Placa do número de série



Fig. 14: Placa do número de série

A placa do número de série está afixada no interior da caixa, do lado direito da traseira do instrumento e apresenta os seguintes dados:

Dados de	Modelo
Identificação	REF: informação de encomenda (número de material e nível de revisão)
	NS: número de série
Dados de endereço	Nome e endereço do fabricante

TECAN



3.3 Dimensões e pesos

	Dimensões	Fluent 480	Fluent 780	Fluent 1080
а	Comprimento total	1150 mm (45,28 in)	1650 mm (64,96 in)	2150 mm (84,65 in)
b	Profundidade da área da base	780 mm (30,71 in)		
с	Profundidade total	ç	923 mm (36,34 in)
d	Altura total no armário	1	1977 mm (77,8 in)

Componente	Fluent 480 Fluent 780		Fluent 1080
Unidade de base	120 kg (264,5 lb) 140 kg (308,6 lb)		190 kg (418,9 lb)
Embalagem	61 kg (135 lb) 83 kg (183 lb)		106 kg (234 lb)
FCA	10,4 kg (22,9 lb)		
RGA	10,2 kg (22,4 lb)		
RGA-Z	10,6 kg (23,4 lb)		
cXP	1,2 kg (2,6 lb)		



Componente	Fluent 480	Fluent 780	Fluent 1080
MCA 96, incluindo o manipulador opcional		19,7 kg (43,43 lb)	

3.4 Fonte de alimentação

AVISO

Sobreaquecimento da fonte de alimentação

A unidade da fonte de alimentação pode ser danificada ou destruída.

- A fonte de alimentação não pode ser tapada.
- É necessário garantir a dissipação do calor da fonte de alimentação.



Não podem ser ligados dispositivos externos à fonte de alimentação. Estes podem resultar na reinicialização ou paragem do Fluent

Tab. 2: Entrada de energia no Fluent

Alimentação	Tensão nominal
Tensão da linha (monofásica)	100–240 V CA
Corrente de entrada	9,8 A (a 100 V) – 4 A (a 240 V)
Frequência	50–60 Hz

Tab. 3: Saída de energia do Fluent

Alimentação	Tensão nominal
Tensão de saída	24–28 V, configuração de fábrica: 25,2 V
Potência contínua	500 W
Potência de pico (limite de tempo)	1500 W durante 3 segundos
Peso	3,8 kg (8,5 lb)

Flutuação máx. da tensão da rede elétrica: ±10% da tensão nominal.

Classificação com vista à segurança elétrica de acordo com as normas EN/CEI:

Tab. 4: Especificações elétricas (segurança)

Categoria de sobretensão	II	IEC 60664-1
Grau de poluição	2	(EN) CEI 61010-1





3.5 Dados e ligações elétricas

A figura apresenta os componentes de sistema de amostra com os dados e as ligações elétricas. Os componentes do instrumento Fluent são apresentados no retângulo. O interruptor de energia do instrumento faz parte da unidade da fonte de alimentação. O fio elétrico é ligado a uma tomada para alimentação de energia da rede.

Todo o tráfego de dados de e para o Fluent passa através da interface USB. O cabo USB é ligado ao PC que controla o instrumento.



3.6 Condições ambientais

▲ CUIDADO

Volumes de pipetagem incorretos

Os resultados da pipetagem podem ser influenciados pelas condições de utilização.

A condensação pode afetar os componentes eletrónicos.

 Se o Fluent for armazenado ou transportado a temperaturas abaixo da temperatura ambiente, irá necessitar de algumas horas para a aclimatação após a instalação.



Condições de utilização O Fluent destina-se apenas à utilização e armazenamento interior.

Temperatura de utilização	15–32 °C (59–90 °F)
Humidade de utilização	30–80% de humidade relativa (sem condensação) a 30 °C (86 °F)
Altitude de utilização	máx. 2000 m acima do nível do mar

Condições de utilização para o manuseio do líquido e a pipetagem:

Temperatura ambiente	20–25 °C (68–77 °F)
Humidade de utilização	30–60% de humidade relativa (sem condensação)
Altitude de utilização	cerca de 500 m acima do nível do mar
Evaporação	Um ambiente com aumento do fluxo de ar (devido ao fluxo laminar, ar condicionado ou ventilação, etc.) aumenta o risco de evaporação, que pode reduzir a precisão da pipetagem, especialmente com volumes baixos ou substâncias voláteis.
	AVISO! Certifique-se de que as condições de validação correspondem às condições da execução.
Temperatura de transporte	-20 a 60 °C (-4 a 140 °F)
Humidade de transporte	20–80% de humidade relativa (sem condensação)
Temperatura de armazenamento	1–60 °C (34–140 °F)

Condições de transporte

Condições de armazenamento CEM



5–80% de humidade relativa (sem condensação) a 30 °C (86 °F) ou nferior

3.7 Emissão e imunidade

Emissão de
ruído< 60 dBA (pressão sonora), medida a uma distância de 1 m do instrumento.
O nível de ruído pode exceder os 78 dB durante breves instantes enquanto o
processo está em execução.

O Fluent está em conformidade com os requisitos de emissão e imunidade descritos na norma IEC 61326-1 e IEC 61326-2-6. No entanto, é necessário avaliar o ambiente eletromagnético antes de utilizar o Fluent. É da responsabilidade do operador garantir a existência de um ambiente eletromagnético compatível ao Fluent para que o Fluent funcione como previsto. O Fluent tem classificação de EQUIPAMENTO DE GRUPO 1 E CLASSE B (CISPR 11).

Este equipamento foi concebido para ser utilizado num AMBIENTE ELETROMAGNÉTICO BÁSICO (IEC 61326-1) e num AMBIENTE DE UNIDADES DE SAÚDE PROFISSIONAIS (IEC 61326-2-6).

É provável que funcione incorretamente se for utilizado num AMBIENTE ELETROMAGNÉTICO INDUSTRIAL (IEC 61326-1) e num AMBIENTE DE UNIDADE DE CUIDADOS DE SAÚDE AO DOMICÍLIO (IEC 61326-2-6). Suspeita-se que o desempenho é afetado por interferências eletromagnéticas; o funcionamento correto pode ser retomado aumentando a distância entre o equipamento e a fonte da interferência.

Não opere o Fluent perto de fontes de radiação eletromagnética forte (por exemplo, fontes de RF intencionais não protegidas), uma vez que podem interferir com a correta operação.

FCC15 Este equipamento foi testado e cumpre os limites previstos para dispositivos digitais de Classe B, de acordo com a parte 15 das Regras da FCC. Estes limites foram concebidos para proporcionar uma proteção razoável contra interferências prejudiciais numa instalação residencial. Este equipamento gera, utiliza e pode irradiar energia de radiofrequência e, se não for instalado e utilizado de acordo com as instruções, pode causar interferências prejudiciais nas comunicações de rádio. No entanto, não há garantia de que não ocorram interferências numa instalação específica. Se este equipamento causar interferências prejudiciais na receção de rádio ou televisão, as quais podem ser determinadas desligando e ligando o equipamento, recomenda-se que o utilizador tente corrigir a interferência através de uma ou mais das seguintes medidas:

- Reorientar ou reposicionar a antena de receção.
- Aumentar a distância entre o equipamento e o recetor.
- Ligar o equipamento a uma tomada num circuito diferente daquele ao qual o recetor está ligado.
- Consulte o revendedor ou um técnico de rádio/TV experiente para obter ajuda.



3.8 Critérios de aceitação da precisão e exatidão da pipetagem

O Fluent Dx só pode ser adquirido com a licença de software Fluent Dx. Não é possível efetuar uma atualização no terreno do instrumento de utilização geral ou do instrumento RUO para Fluent Dx.

Se uma opção ou dispositivo para utilização apenas em investigação (RUO) estiver integrado com o Fluent, a utilização prevista muda para apenas para utilização em investigação. Não deve ser utilizado em procedimentos de diagnóstico.

6

Ĭ

1

Uma integração no local do MCA 384 no Fluent Dx irá alterar a utilização prevista para "Utilização geral" e o manual de utilização (ID doc. 399706) será aplicado.

Tab. 5: Critérios de aceitação da precisão e exatidão da pipetagem Liquid Flexible Channel Arm (Liquid FCA) para utilização com a deteção capacitiva do nível do líquido (cLLD)

Classe de líquidos cLLD do Fluent Control	Liquid FCA								
	Ponta de perfuração , seringa de 5000 ul, tubo de volume padrão	Ponta de perfuração , seringa de 1250 ul, tubo de volume padrão	Ponta fixa padrão, seringa de 1250 ul, tubo de volume padrão	Ponta de volume reduzido, seringa de 250 ul, tubo de volume reduzido	Te-PS, seringa de 250 ul, tubo de volume reduzido				
DMSO, contacto húmido simples			Exemplo	Exemplo	Exemplo				
DMSO, livre, diversos			12 x 50 ul CV≤1,0% AC±2,0%	12 x 5 ul CV≤3,5% AC±5,0%	Exemplo				
DMSO, livre, simples			5 ul CV≤3,0% AC±6,0%	1 ul CV≤4,0% AC±5,0%	1 ul CV≤5,0% AC±8,0%				
			10 ul CV≤2,0% AC±2,0%	2 ul CV≤3,5% AC±4,0%	2 ul CV≤3,0% AC±6,0%				
			1000 ul CV≤0,25% AC±0,75%	10 ul CV≤2,0% AC±2,0%	10 ul CV≤0,8% AC±1,0%				



Tab.	6: Critérios c	le aceitação d	a precisá	ăo e exati	dão da pi	petagem	Liquid Flex	ible Cha	annel
Arm	(Liquid FCA)	para utilizaçã	o com a	deteção d	capacitiva	do nível	do líquido ((cLLD)	

Classe de líquidos cLLD do Fluent Control	Liquid FCA								
	DiTi 1000, abertur a larga, seringa de 1250 ul	DiTi 1000, seringa de 1250 ul, tubo de volume padrão	DiTi 350, seringa de 1250 ul, tubo de volume padrão	DiTi 200, seringa de 1250 ul, tubo de volume padrão	DiTi 50, seringa de 1250 ul, tubo de volume padrão	DiTi 10, seringa de 1250 ul, tubo de volume padrão			
DMSO, contacto húmido simples		Exemplo		Exemplo	1000 ul CV≤8,0% AC±10,0%				
DMSO, livre, diversos		12 x 50 ul CV≤2,5% AC±3,0%	Exemplo	12 x 10 ul CV≤6,0% AC±4,0%	6 x 5 ul CV≤5,0% AC±5,0%				
DMSO, livre, simples		10 ul 10 u CV≤2,5% CV AC±3,0% AC	10 ul CV≤2,0% AC±2,0%	10 ul CV≤2,0% AC±2,0%	1 ul CV≤8,0% AC±10,0%	1 ul CV≤8,0% AC±10,0%			
					2 ul CV≤3,0% AC±5,0%				
		100 ul CV≤0,5% AC±1,5%	100 ul CV≤0,4% AC±1,5%	200 ul CV≤0,4% AC±1,0%	10 ul CV≤1,5% AC±2,0%	10 ul CV≤2,0% AC±2,0%			
					50 ul CV≤0,75% AC±2,0%				

Tab. 7: Critérios de aceitação da precisão e exatidão da pipetagem Liquid Flexible Channel Arm (Liquid FCA) para utilização com a deteção capacitiva do nível do líquido (cLLD)

Classe de líquidos cLLD do Fluent Control	Liquid FCA								
	Ponta de perfuração , seringa de 5000 ul, tubo de volume padrão	Ponta de perfuração , seringa de 1250 ul, tubo de volume padrão	Ponta fixa padrão, seringa de 1250 ul, tubo de volume padrão	Ponta de volume reduzido, seringa de 250 ul, tubo de volume reduzido	Te-PS, seringa de 250 ul, tubo de volume reduzido				
Etanol, livre, diversos			12 x 50 ul³ CV≤2,5% AC±3,0%						



Classe de líquidos cLLD do Fluent Control	Liquid FCA							
	Ponta de perfuração , seringa de 5000 ul, tubo de volume padrão	Ponta de perfuração , seringa de 1250 ul, tubo de volume padrão	Ponta fixa padrão, seringa de 1250 ul, tubo de volume padrão	Ponta de volume reduzido, seringa de 250 ul, tubo de volume reduzido	Te-PS, seringa de 250 ul, tubo de volume reduzido			
Etanol, livre, simples			50 ul³ CV≤1,0% AC±3,0%					
Soro, livre, diversos			Exemplo					
Soro, livre, simples			5 ul CV≤5,0% AC±8,0%					
Sangue total com perfuração simples ⁵		200 ul ⁶ CV≤2,0% AC±3,0%						
Sangue total com diversas perfuraçõe s	4 x 1000 ul CV≤1,5% AC±2,0%							

Tab. 8: Critérios de aceitação da precisão e exatidão da pipetagem Liquid Flexible Channel Arm (Liquid FCA) para utilização com a deteção capacitiva do nível do líquido (cLLD)

Classe de líquidos cLLD do Fluent Control	Liquid FCA								
	DiTi 1000, abertura larga, seringa de 1250 ul	DiTi 1000, seringa de 1250 ul, tubo de volume padrão	DiTi 350, seringa de 1250 ul, tubo de volume padrão	DiTi 200, seringa de 1250 ul, tubo de volume padrão	DiTi 50, seringa de 1250 ul, tubo de volume padrão	DiTi 10, seringa de 1250 ul, tubo de volume padrão			
Etanol, livre, diversos				6 x 10 ul³ CV≤6,0% AC±4,0%					
Etanol, livre, simples		Exemplo		Exemplo	30 ul ³ CV≤0,75% AC±2,0%				



Classe	Liquid FCA								
de líquidos cLLD do Fluent Control	DiTi 1000, abertura larga, seringa de 1250 ul	DiTi 1000, seringa de 1250 ul, tubo de volume padrão	DiTi 350, seringa de 1250 ul, tubo de volume padrão	DiTi 200, seringa de 1250 ul, tubo de volume padrão	DiTi 50, seringa de 1250 ul, tubo de volume padrão	DiTi 10, seringa de 1250 ul, tubo de volume padrão			
Soro, livre, diversos	12 x 50 ul CV≤5,5% AC±5,0%	Exemplo	Exemplo	6 x 25 ul CV≤3,5% AC±5,0%	Exemplo				
	6 x 100 ul CV≤3,5% AC±5,0%								
	4 x 200 ul CV≤3,5% AC±5,0%								
Soro, livre, simples	20 ul CV≤5,5% AC±8,0%	Exemplo	Exemplo	Exemplo	10 ul CV≤2,0% AC±5,0%	Exemplo			
	100 ul CV≤1,5% AC±2,0%								
	500 ul CV≤0,5% AC±1,0%								
	900 ul CV≤0,5% AC±1,0%								
Sangue total com perfuraç ão simples									
Sangue total com diversas perfuraç ões									



Tab. 9: Critérios de aceitação da precisão e exatidão da pipetagem Liquid Flexible Channel Arm (Liquid FCA) para utilização com a deteção capacitiva do nível do líquido (cLLD)

Classe de líquidos cLLD do Fluent Control	Liquid FCA								
	Ponta de perfuração , seringa de 5000 ul, tubo de volume padrão	Ponta de perfuração , seringa de 1250 ul, tubo de volume padrão	Ponta fixa padrão, seringa de 1250 ul, tubo de volume padrão	Ponta de volume reduzido, seringa de 250 ul, tubo de volume reduzido	Te-PS, seringa de 250 ul, tubo de volume reduzido				
Água, perfuração simples		200 ul ⁶ CV≤1,5% AC±3,0%							
Água, contacto húmido simples	Exemplo	Exemplo	Exemplo	Exemplo	0,2 ul CV ≤10,0% AC±15,0% (valor não incluído no LC)				
Água, livre, diversos			12 x 50 ul CV≤2,0% AC±2,0%	12 x 5 ul CV≤2,5% AC±5,0%	Exemplo				
Água, livre, simples			5 ul CV≤3,0% AC±3,0%	0,5 ul CV≤6,0% AC±10,0%	0,5 ul CV≤6,0% AC±10,0%				
			10 ul CV≤1,75% AC±2,5%	1 ul CV≤3,5% AC±8,0%	1 ul CV≤4,0% AC±6,0%				
			1000 ul CV≤0,3% AC±0,75%	10 ul CV≤1,0% AC±2,0%	10 ul CV≤0,8% AC±1,0%				



Tab. 10: Critérios de aceitação da precisão e exatidão da pipetagem Liquid Flexible Channel Arm (Liquid FCA) para utilização com a deteção capacitiva do nível do líquido (cLLD)

Classe	Liquid FCA								
de líquido s cLLD do Fluent Contro I	DiTi 1000, abertura larga, seringa de 1250 ul	DiTi 1000, seringa de 1250 ul, tubo de volume padrão	DiTi 350, seringa de 1250 ul, tubo de volume padrão	DiTi 200, seringa de 1250 ul, tubo de volume padrão	DiTi 50, seringa de 1250 ul, tubo de volume padrão	DiTi 10, seringa de 1250 ul, tubo de volume padrão			
Água, perfura ção simple s									
Água, contact o húmido simple s		Exemplo		Exemplo	Exemplo	Exemplo			
Água, livre, diverso s	6 x 100 ul CV≤6,5% AC±5,0%	12 x 50 ul CV≤2,0% AC±2,0%	12 x 50 ul CV≤2,0% AC±2,0%	12 x 10 ul CV≤3,5% AC±2,0%	12 x 10 ul CV≤3,5% AC±1,5%	6 x 5 ul CV≤4,0% AC±3,0%			
	4 x 200 ul CV≤3,5% AC±5,0%								
Água, livre, simple	20 ul CV≤2,5% AC±5,0%	10 ul CV≤2,0% AC±3,0%	10 ul CV≤2,0% AC±2,0%	10 ul CV≤1,8% AC±2,5%	1 ul CV≤6,0% AC±8,0%	0,5 ul CV≤8,0% AC±10,0%			
S	100 ul CV≤1,0% AC±2,0%	100 ul CV≤0,3% AC±0,5%			2 ul CV≤3,0% AC±4,0%	1 ul CV≤4,0% AC±5,0%			
	500 ul CV≤0,5% AC±2,0%	1000 ul CV≤0,2% AC±0,5%	100 ul CV≤0,3% AC±1,0%	200 ul CV≤0,2% AC±0,75%	10 ul CV≤1,0% AC±2,0%	10 ul CV≤1,0% AC±1,0%			
	900 ul CV≤0,5% AC±2,0%				50 ul CV≤0,5% AC±1,5%				



Tab. 11: Critérios de aceitação da precisão e exatidão da pipetagem Air Flexible Channel Arm MultiSense (Air FCA MS) para utilização com a deteção capacitiva do nível do líquido (cLLD)

Classe de líquidos cLLD do Fluent Control	Air FCA							
	DiTi 1000 ul, abertura larga	DiTi 1000 ul	DiTi 1000 ul transparente	DiTi 350 ul				
DMSO, contacto húmido simples								
DMSO, livre, diversos		Exemplo		Exemplo				
DMSO, livre, simples		Exemplo		Exemplo				

Tab. 12: Critérios de aceitação da precisão e exatidão da pipetagem Air Flexible Channel Arm MultiSense (Air FCA MS) para utilização com a deteção capacitiva do nível do líquido (cLLD)

Classe de líquidos cLLD do Fluent Control	Air FCA							
	DiTi 200 ul	DiTi 200 ul transparen te	DiTi 50 ul	DiTi 50 ul transparen te	DiTi 10 ul			
DMSO, contacto húmido simples			1 ul CV≤8,0% AC±10,0%					
DMSO, livre, diversos	Exemplo		4 x 10 ul CV≤5,0% AC±5,0%					
DMSO, livre, simples	10 ul CV≤1,5% AC±3,0%		1 ul CV≤6,0% AC±5,0%		Exemplo			



Tab. 13: Critérios de aceitação da precisão e exatidão da pipetagem Air Flexible Channel Arm MultiSense (Air FCA MS) para utilização com a deteção capacitiva do nível do líquido (cLLD)

Classe de	Air FCA			
cLLD do Fluent Control	DiTi 1000 ul, abertura larga	DiTi 1000 ul	DiTi 1000 ul transparente	DiTi 350 ul
Etanol, livre, diversos		Exemplo		Exemplo
Etanol, livre, simples		Exemplo 1 C A 5 C A	100 ul CV≤1,5% AC±2,5%	Exemplo
			500 ul CV≤1,0% AC±2,5%	
MasterMix, livre, diversos		Exemplo		Exemplo
MasterMix, livre, simples		Exemplo		Exemplo
Soro, livre, diversos	12 x 25 ul CV≤8,0% AC±5,0%	6 x 100 ul CV≤2,0% AC±1,0%		Exemplo
	12 x 50 ul CV≤5,0% AC±5,0%			
	6 x 100 ul CV≤3,0% AC±5,0%			
	4 x 200 ul CV≤3,0% AC±5,0%			
Soro, livre, simples	20 ul CV≤5,0% AC±5,0%	Exemplo	10 ul CV≤6% AC±5,0%	100 ul CV≤0,3% AC±0,5%
	100 ul CV≤1,0% AC±2,0%	100 ul CV≤1,0% AC±2,0%		
	500 ul CV≤0,5% AC±1,0%	ıl),5% I,0%	500 ul CV≤0,5% AC±2,0%	
900 ul CV≤0,5% AC±1,0%				



Tab. 14: Critérios de aceitação da precisão e exatidão da pipetagem Air Flexible Channel Arm MultiSense (Air FCA MS) para utilização com a deteção capacitiva do nível do líquido (cLLD)

Classe de	Air FCA				
cLLD do Fluent Control	DiTi 200 ul	DiTi 200 ul transparen te	DiTi 50 ul	DiTi 50 ul transparen te	DiTi 10 ul
Etanol, livre, diversos	6 x 10 ul CV≤3,0% AC±4,0%		Exemplo		
Etanol, livre, simples	Exemplo	3 ul CV≤6,0% AC±12,0%	40 ul CV≤1,0% AC±1,0%	1 ul CV≤8,0% AC±15,0%	Exemplo
		5 ul CV≤6,0% AC±12,0%		10 ul CV≤3,0% AC±4,0%	
	_	10 ul CV≤3,5% AC±5,0%		40 ul (com filtro) CV≤1,0% AC±2,0%	
		100 ul CV≤1,0% AC±4,0%		50 ul (sem filtro) CV≤1,0% AC±2,0%	
MasterMix, livre, diversos	Exemplo		6 x 5 ul CV≤5,0% AC±5,0%		
		4 x 10 ul CV≤3,0% AC±5,0%			
MasterMix, livre, simples	Exemplo		5 ul CV≤3,0% AC±5,0%		Exemplo
Soro, livre, diversos	6 x 5 ul CV≤8,0% AC±7,5%		6 x 5 ul CV≤8,0% AC±5,0%		
	6 x 25 ul CV≤3,0% AC±3,0%				



Classe de	Air FCA				
cLLD do Fluent Control	DiTi 200 ul	DiTi 200 ul transparen te	DiTi 50 ul	DiTi 50 ul transparen te	DiTi 10 ul
Soro, livre, simples	100 ul CV≤0,3% AC±0,5%	5 ul CV≤3% AC±8,0%	5 ul CV≤3,5% AC±5,0%		Exemplo
		10 ul CV≤2,0% AC±6,0%	10 ul CV≤1,0% AC±1,0%	-	
	100 ul CV≤1,0% AC±2,0%				

Tab. 15: Critérios de aceitação da precisão e exatidão da pipetagem Air Flexible Channel Arm MultiSense (Air FCA MS) para utilização com a deteção capacitiva do nível do líquido (cLLD)

Classe de	Air FCA				
cLLD do Fluent Control	DiTi 1000 ul, abertura larga	DiTi 1000 ul	DiTi 1000 ul transparente	DiTi 350 ul	
Água, livre, diversos	6 x 100 ul CV≤6,0% AC±5,0%	12 x 50 ul CV≤2,0% AC±1,0%		12 x 10 ul CV≤4,0% AC±2,0%	
	4 x 200 ul CV≤3,0% AC±5,0%	6 x 100 ul CV≤1,0% AC±1,5%		6 x 20 ul CV≤1,5% AC±2,5%	
Água, livre, simples	20 ul CV≤2,0% AC±5,0%	10 ul CV≤1,2% AC±2,0%	10 ul CV≤3% AC±8,0%	10 ul CV≤0,6% AC±1,5%	
	100 ul CV≤0,5% AC±2,0%		100 ul CV≤0,75% AC±2,0%		
	500 ul CV≤0,5% AC±2,0%	100 ul CV≤0,3% AC±0,5%	500 ul CV≤0,5% AC±1,0%	200 ul CV≤0,2% AC±0,5%	
	900 ul CV≤0,5% AC±2,0%				



Tab. 16: Critérios de aceitação da precisão e exatidão da pipetagem Air Flexible Channel Arm MultiSense (Air FCA MS) para utilização com a deteção capacitiva do nível do líquido (cLLD)

Classe de	Air FCA					
cLLD do Fluent Control	DiTi 200 ul	DiTi 200 ul transparen te	DiTi 50 ul	DiTi 50 ul transparen te	DiTi 10 ul	
Água, livre, diversos	12 x 10 ul CV≤4,0% AC±2,0%		6 x 5 ul CV≤6,0% AC±5,0%			
Água, livre, simples 10 ul CV≤0,6% AC±1,5%	10 ul CV≤0,6% AC±1,5%	5 ul CV≤4% AC±8,0%	1 ul CV≤4,0% AC±8,0%	1 ul CV≤10,0% AC±20,0%	0,5 ul CV≤5,0% AC±9,5%	
		10 ul CV≤3,0% AC±4,0%	5 ul CV≤0,8% AC±1,5%	10 ul CV≤1,5% AC±2,0%		
200 ul CV≤0, AC±0,	200 ul 100 ul CV≤0,2% CV≤0,75% AC±0,5% AC±1,0%	10 ul CV≤0,5% AC±1,0%	40 ul (com filtro) CV≤0,5% AC±1,0%	10 ul CV≤ 0,8% AC±2,0%		
			50 ul CV≤0,3% AC±0,5%	50 ul (sem filtro) CV≤0,5% AC±1,0%		

Tab. 17: Critérios de aceitação da precisão e exatidão da pipetagem Air Flexible Channel Arm MultiSense (Air FCA MS) para utilização com a deteção de pressão do nível do líquido (pLLD)

Classe de	Air FCA				
pLLD do FluentContro	DiTi1000F	DiTi200F	DITI50F ¹	DiTi10F	
Hexano, livre, simples	100 ul CV≤1,0% AC±1,5%	10 ul CV≤3,5% AC±2%	10 ul CV≤3% AC±1,5%	Exemplo	
	500 ul CV≤0,6% AC±1,0%	100 ul CV≤1,5% AC±1%	45 ul CV≤1,5% AC±1,0%		





Classe de	Air FCA				
pLLD do FluentContro I	DiTi1000F	DiTi200F	DiTi50F ¹	DITI10F	
Água, livre, simples²	10 ul CV≤2% AC±2%	10 ul CV≤2% AC±2%	5 ul CV≤1,5% AC±1,5%	2 ul CV≤5% AC±9,5%	
	100 ul CV≤0,5% AC±1%	195 ul CV≤0,5% AC±1%	10 ul CV≤2% AC±2,5%	10 ul CV≤ 2% AC±3%	

Tab. 18: Critérios de aceitação da precisão e exatidão da pipetagem Multi Channel Arm 96 (MCA 96) para utilização com a deteção capacitiva do nível do líquido (cLLD)

Classe de líquidos	MCA 96		
CLLD do FluentControl	DiTi 1000 F	DiTi 10F	
DMSO, livre, simples	1000 ul CV≤1,5%	1 ul CV≤9%	
Água, livre, simples	1000 ul CV≤1,5%	1 ul CV≤4%	

Chave de especificação de desempenho (exemplo)

Volume mais baixo testado	10 ul
Especificação de CV	CV≤2,0%
Especificação de exatidão	AC±3%

Definições

• CV:

medição da precisão de distribuição que indica a extensão da variabilidade em relação à média das medições.

• AC:

exatidão da distribuição que indica a proximidade da medição com o valor pretendido.

Condições de medição

Os valores foram medidos de acordo com o seguinte protocolo.

- Medições obtidas em instrumentos Fluent padrão operados de acordo com as condições ambientais especificadas para o manuseamento de líquidos e mantidos de acordo com as instruções de cuidados do sistema.
- Mínimo de 96 medições
- A CV especificada e a exatidão média são os valores máximos obtidos no geral e por canal.





Apenas as pontas descartáveis Tecan garantem o desempenho especificado para os instrumentos de pipetagem Tecan.

Pode ser necessária calibração para volumes de 5 µl ou inferiores.

Os valores de pipetagem de DMSO são obtidos com água como líquido do sistema.

É utilizada uma formulação MasterMix com 50% de glicerol para o Air FCA.

Notas

- ¹ Especificado para o DiTi de 50 ul (número da peça 30200712) (específico do fornecedor)
- ² Classe de líquido pLLD de água, livre, simples, suportado a partir da versão 3.7 do FluentControl
- ³ Obtido com o comando de pré-recolha/obtenção do bloco de pontas para expelir a solução de lavagem
- ⁴ Possibilidade de 0,2 ul com personalização avançada de parâmetros (contactar linha especializada para obter informações)
- ⁵ Com volume excessivo
- ⁶ Especificações de desempenho para distribuição livre em material de laboratório não perfurável
- Exem Modelo de classe de líquido fornecido no FluentControl sem plo especificação de desempenho

Célul Nenhum modelo de classe de líquido fornecido no FluentControl a em

branc o

ul Forma alternativa para microlitro, utilizada em vez de µl

Média



$$\overline{x} = \frac{1}{n} \sum_{i=1}^{n} x_i$$

Desvio-padrão

$$s = \sqrt{\frac{1}{n-1} \sum_{i=1}^{n} (x_i - \overline{x})^2}$$

Coeficiente de variação

$$CV = \frac{s}{x} * 100\%$$

Exatidão





- x Média
- n Número de pontos de dados
- Σ Símbolo de soma
- I Indicador a começar em 1
- x_i i valor do conjunto de dados
- s Desvio-padrão
- CV Coeficiente de variação

Teste de contaminação cruzada

Método de tabuleiro de xadrez para avaliação de contaminação cruzada

Critérios de aceitação:

os valores de Ct dos corantes fluorescentes FAM e VIC no controlo sem modelo (NTC) ou nos poços de controlo negativo têm de ser N/A ou >37 para todos os processamentos. Além disso, os valores de Ct de FAM e VIC nos poços de controlo positivo têm de ser válidos (≤37). O valor de Ct indica o número de ciclo em que o sinal fluorescente de uma amostra excede um limite, indicando a deteção do ácido nucleico alvo.

Resultados:

abaixo é fornecido um exemplo de resultados de qPCR. O primeiro gráfico mostra as curvas de amplificação esperadas dos poços de controlo positivo, com valores de Ct <37. O segundo gráfico confirma a ausência de amplificação nos poços de controlo sem modelo, com valores de Ct registados como N/A.







4 Descrição do funcionamento

Este capítulo explica o funcionamento básico do Fluent, ilustra a estrutura e fornece uma descrição funcional dos elementos.

4.1 Vista geral

O Fluent é utilizado para tarefas de pipetagem com braços robóticos. Os braços robóticos podem aspirar de e distribuir para vários recipientes, tais como tubos de ensaio ou microplacas.

O Fluent está disponível em três tamanhos diferentes:

- Fluent 480
- Fluent 780
- Fluent 1080



Fig. 16: Vista geral do instrumento (o instrumento pode diferir da ilustração)

A Plataforma

- Braços robóticos
- **C** Opções e dispositivos



Também está disponível um conjunto de proteção antissísmica para áreas de alto risco sísmico.

В

Para obter mais informações, consulte a secção "Apoio ao cliente" [> 213].

4.2 Mesa de trabalho

Segmentos A mesa de trabalho do Fluent, que é a zona de amostras do instrumento, é composta por segmentos. Os segmentos da mesa de trabalho são componentes da mesa de trabalho intercambiáveis que podem ter várias dimensões e características. CUIDADO! Não opere o sistema sem os segmentos da mesa de trabalho.
Número da A largura do segmento é expressa em números de grelha. A grelha tem 25 mm de

nero da
grelhaA largura do segmento é expressa em números de grelha. A grelha tem 25 mm de
largura e corresponde à distância entre os pinos de posicionamento de um
segmento.





Os números da grelha também são utilizados para expressar a localização dos segmentos ou calhas da mesa de trabalho.

4.2.1 Carregadores



Fig. 17: Mesa de trabalho do Fluent



Os carregadores são componentes da mesa de trabalho concebidos para conter o material laboratorial ou os consumíveis na mesa de trabalho.

As calhas são carregadores que deslizam para dentro e para fora dos segmentos da grade e normalmente contêm os tubos de ensaio ou cartuchos de reagente.

Os segmentos são elementos estáticos, bloqueados na mesa de trabalho. Alguns segmentos têm recetáculos (segmentos de recetáculo) que contêm material laboratorial, como as microplacas ou as placas deep well, ou consumíveis como as caixas DiTi. Alguns segmentos têm pinos de grelha (segmentos da grade) para carregar e descarregas as calhas.





4.2.2 Tabuleiros de recolha da mesa de trabalho

Fig. 18: Tabuleiro de recolha

Os tabuleiros de recolha, que são colocados por baixo dos segmentos dinâmicos da plataforma, recolhem os derrames de líquido que podem ocorrer na área de carregamento manual da plataforma. O sistema deve ser operado com o maior número possível de tabuleiros de recolha instalados sob a plataforma para recolher todos os derrames de líquidos. **CUIDADO! Não operar o sistema sem os tabuleiros de recolha e os segmentos da plataforma.**

Apenas a versão armário pode ter recortes nos tabuleiros de recolha para ferramentas e instrumentos.

As áreas elevadas em cada tabuleiro de recolha foram concebidas para permitir recortes sem afetar o volume de captação. Estes recortes acomodam calhas para resíduos ao longo da plataforma ou outras integrações de dispositivos ao utilizar a versão montada do armário. Para os pontos de acesso ao longo da plataforma, os tabuleiros podem ser colocados numa prateleira no armário, por baixo dos dispositivos integrados.



Fig. 19: Tabuleiros de recolha por baixo dos segmentos da mesa de trabalho

Os tabuleiros de recolha não estarão presentes nos casos em que o RGA necessitar de aceder a um dispositivo por baixo da plataforma. Está incluído com o instrumento um conjunto de tabuleiros de recolha. Os tabuleiros de recolha podem ser lavados ou substituídos, conforme necessário. Consulte a secção Final do dia.



4.2.3 Colocação no segmento



Fig. 20: Pinos de bloqueio e pinos de posicionamento

A Pinos de bloqueio B Pinos de posicionamento

O Fluent utiliza pinos para posicionar corretamente as calhas, os adaptadores ou as opções num segmento. As calhas destinam-se a deslizar nos pinos. O seu posicionamento pode ser verificado pela leitura do número da grelha, na parte da frente do instrumento. Os pinos de bloqueio fixam as calhas na posição.





4.2.4 Posição do segmento

Fig. 21: Posições laterais e posições de grelha

A Posições laterais

As posições laterais (AB, YZ) podem ser utilizadas para colocar o material laboratorial manuseado pelo RGA.

В

Posições de grelha



Com o FCA ou MCA, não é possível realizar a pipetagem em posições laterais.

Os braços de pipetagem podem aceder às posições de grelha numeradas (1–n). Com múltiplas configurações de braços, no entanto, nem todas as grelhas numéricas estão acessíveis a todos os braços de pipetagem. Podem aplicar-se restrições, dependendo da configuração do braço do instrumento.

4.3 Braços robóticos

O Fluent pode ser equipado com diversos braços robóticos:

- Flexible Channel Arm (FCA)
- Multiple Channel Arm (MCA 96)
- Robotic Gripper Arm (RGA)

Os braços robóticos podem ser equipados com vários acessórios para braços.



4.3.1 Flexible Channel Arm (FCA)



Fig. 22: Flexible Channel Arm



Se não forem tolerados resíduos de transição, recomendamos vivamente a utilização de pontas descartáveis com filtro.

O FCA (A) está equipado com pontas de pipetagem e pode controlar o manuseio do líquido para até 8 canais separados.

O FCA configurado com adaptadores DiTi possui um manipulador FCA opcional que permite determinados movimentos de material laboratorial—consulte "Manipulador FCA" [> 75].

4.3.1.1 FCA com sistema de líquido (Liquid FCA)

O FCA com um sistema de deslocamento de líquido está cheio com líquido do sistema que é fornecido pelas bombas de seringa. É utilizado para pipetar líquidos com diferentes âmbitos de volume, dependendo das pontas e da seringa utilizadas. O Liquid FCA pode ser configurado com uma seleção de pontas fixas laváveis ou com adaptadores de pontas descartáveis.



A Tecan recomenda a utilização de água desionizada como líquido do sistema.

4.3.1.2 FCA com Sistema de ar (Air FCA)

O FCA com um sistema de deslocamento de ar é utilizado para pipetar líquidos, através do deslocamento de um êmbolo no interior do canal de pipetagem. O Air FCA está configurado com adaptadores de pontas descartáveis.

4.3.2 Multiple Channel Arm 96 (MCA 96)

O MCA 96 é um braço robótico com uma cabeça de pipetagem com múltiplos canais. Os 96 canais da cabeça de pipetagem aspiram e distribuem em simultâneo. O braço MCA 96 tem as seguintes caraterísticas principais:





Fig. 23: Cabeça/Manipulador do MCA 96

- · Compatível com o portefólio de DiTi condutoras do FCA até 1000 µl, inclusive
- Intervalo de pipetagem de 1 µl a 1000 µl
- Efetua uma deteção (capacitiva) do nível de líquido (cLLD)
- Efetua uma recolha parcial de DiTi e uma pipetagem de compensação (por exemplo, para diluição) ao utilizar produtos de pontas com tabuleiro combinado Tecan
- Manipulador opcional para tarefas simples de manuseamento de material laboratorial



A cabeça do MCA 96 funciona com pontas descartáveis condutoras no MCA 96 para compatibilidade com a cLLD.

A cLLD no MCA 96 apenas está verificada para ser utilizada em recipientes.

A primeira ponta a entrar num líquido aciona o sinal da cLLD. Alturas de enchimento desiguais (por exemplo, em poços de uma microplaca enchidos manualmente) podem levar a um processamento não intencional e a resultados erróneos.





Ao utilizar o MCA 96, utilizar apenas material laboratorial e suportes compatíveis com cLLD. Utilizar tabuleiros compatíveis para a recolha parcial de DiTi: Tabuleiro SLAS Tecan. Para obter mais informações, consultar o manual de referência (ref.^a [4]).

O manipulador opcional do MCA 96, atualizável em campo, cumpre a finalidade de tarefas simples de manuseamento de material laboratorial.

O manipulador pode ser utilizado, por exemplo, para as seguintes tarefas:

- Mover as microplacas e as DWP no modo horizontal para as posições de destino na mesa de trabalho
- Trabalhar com tabuleiros de DiTi empilhados, ao descartar camadas vazias para o lixo
- Manusear tampas de microplacas

O manipulador do MCA foi concebido como um submódulo na cabeça do MCA, pelo que é uma alternativa compacta e económica a um braço manipulador robótico (RGA) dedicado. O manipulador do MCA 96 não é capaz de efetuar tarefas de manuseamento de material laboratorial mais complexas, tais como:

- Acesso a hotéis
- Manuseamento do material laboratorial na orientação vertical
- Tarefas que incluem rotações

Estas tarefas complexas ainda requerem a adição de um braço manipulador robótico (RGA) dedicado ao sistema. O manipulador do MCA 96 é composto pela mesma estação de troca de dedos que o manipulador RGA.

4.3.3 Robotic Gripper Arm (RGA)



Fig. 24: Robotic Gripper Arm

O RGA (A) é um braço robótico com uma cabeça do manipulador e dedos do manipulador. O RGA transporta as microplacas e outro material laboratorial entre as posições da mesa de trabalho, os dispositivos periféricos e o armazenamento do material laboratorial:

Um Robotic Gripper Arm com altura standard (RGA Z standard) pode aceder a objetos localizados na mesa de trabalho ou na plataforma inferior.



Um Robotic Gripper Arm alto (RGA Z longo) pode aceder a objetos localizados na mesa de trabalho, na mesa de trabalho inferior e abaixo.

4.3.3.1 Cabeça do manipulador robótico

O RGA pode ser equipado com duas opções diferentes de cabeça do manipulador robótico.

A cabeça do manipulador normal oferece uma seleção de dedos do manipulador que são permutáveis manualmente.

O Sistema de troca de dedos (FES) oferece a troca automática de dedos com uma seleção de conjuntos de dedos do manipulador. Os conjuntos de dedos são montados numa estação de encaixe montada num segmento de recetáculo standard. Os conjuntos de dedos são automaticamente recolhidos e colocados pelo braço robótico. O tipo de dedo e a troca de dedos são monitorizados. Num único método, é possível utilizar qualquer dedo ou todos os dedos.

4.3.4 Acessórios do braço



D

Fig. 25: Acessórios do braço

- A Pontas fixas
- **C** Dedos do manipulador
- B Pontas descartáveis
 - Leitor de código de barras para o RGA

4.3.4.1 Pontas fixas

Se não forem tolerados resíduos de transição, recomendamos vivamente a utilização de pontas descartáveis com filtro.

Estão disponíveis pontas reutilizáveis laváveis para a aspiração e distribuição para o FCA e o MCA.

4.3.4.2 Pontas descartáveis

As pontas são disponibilizadas em tabuleiros ou caixas (individuais ou agrupadas) dependendo do tipo. As pontas são eliminadas ou colocadas novamente no rack após a aspiração. As pontas são eliminadas, com o Sistema de ejeção de pontas descartáveis, numa rampa de resíduos montada num segmento da mesa de trabalho.



4.3.4.3 Dedos do manipulador

Estão disponíveis diferentes tipos de dedos do manipulador para a cabeça do manipulador RGA normal e para o manipulador do MCA 96, ambos com o sistema de troca de dedos (FES).

Dedos doOs dedos do manipulador excêntricos transportam objetos baseados emmanipuladormicroplacas dentro e além da área de pipetagem. Agarram objetos das placas aexcêntricospartir dos lados. Estão disponíveis duas variantes:

- Dedos com comprimento padrão para carregar as microplacas para hotéis e dispositivos.
- Dedos excêntricos longos para carregar dispositivos mais profundos, como a incubadora monitorizada de placa de células com 4 ranhuras.

Dedos do manipulador excêntricos do MCA 96

> Dedos do manipulador

> > cêntricos

Os dedos do manipulador excêntricos do MCA 96 transportam objetos à base de microplacas dentro da área de pipetagem. Este tipo de dedo consegue agarrar objetos da placa por baixo da cabeça do manipulador do MCA 96. (Este tipo de dedo NÃO é compatível com o manipulador RGA.)

Os dedos do manipulador cêntricos transportam objetos baseados em microplacas dentro e abaixo da área de pipetagem. Agarram objetos das placas a partir de cima. Aplicável apenas para a cabeça do manipulador RGA normal.

Dedos de tubo Os dedos de tubo transportam objetos baseados em tubos dentro e abaixo da área de pipetagem. Aplicável apenas para a cabeça do manipulador RGA normal.

4.3.4.4 Leitor de código de barras

O RGA pode ser equipado com um leitor horizontal para os códigos de barras das microplacas e caixas DiTi.



As instruções de segurança da classe laser devem ser lidas atentamente e seguidas.

Consulte também o manual disponibilizado pelo fabricante do leitor de códigos de barras.

4.4 Sistema de líquido (Liquid FCA)



Fig. 26: Sistema de líquido (Liquid FCA)

O sistema de líquido destina-se à lavagem eficiente, no interior e exterior, das pontas de pipetagem fixas.


4.5 Opções e dispositivos

	Nesta secção, as opções e os dispositivos para utilização apenas em investigação encontram-se marcados com um asterisco (*).			
	Para obter mais informações, consulte a secção "Utilização prevista" [> 8].			
Opções	Hotel (dispositivo de armazenamento de placas)			
passivas	Armário			
	Cobertura contra pó			
	Manipulador FCA			
pções ativas	Filtro HEPA			
	Fluent Stacker			
	• MIO2			
	• Te-Shake			
	• Te-VacS			
	Tapete rolante Eluent			
	Resolvex i300			
	Pontas de perfuração e rotador de tubos			
	Leitor FRIDA			
	Para obter mais informações, consulte a secção "Documentos de referência" [▶ 9].			
	 Dispositivos de lavagem baseados no HydroControl 			
	Balanças baseadas na norma MT-SICS nível 1			
	 Dispositivos compatíveis com a SiLA* 			
	Selador Agilent*			
	ODTC Inheco			
	Aquecimento e arrefecimento Inheco, utilizando o controlador MTC/STC			
	• Cytomat 10*, 20*, 200* e 6000*			
Leitores de	Leitor de códigos de barras de tubos Fluent ID			
códigos de barras	Leitores de códigos de barras da série Keynence BL-1300			
Leitores	Leitores Tecan controlados pelo Magellan			
	Spark e SparkControl Magellan*			
	 Leitor plano Ziath 2D* 			

Fluent® Manual de utilização, 403096, pt, V1.0





4.5.1 Leitor de códigos de barras de tubos Fluent ID

Fig. 27: Fluent ID

- A Caixa do leitor
- **C** Área de carregamento
- B Leitor de código de barras a laserD LED

E Refletor

O Fluent ID é um módulo opcional que pode ser incorporado para ler as etiquetas de códigos de barras dos tubos à medida que as calhas dos tubos são carregadas na mesa de trabalho. Cada módulo do Fluent ID inclui seis posições de grelha dedicadas para carregar e ler as etiquetas de códigos de barras de até seis calhas. É utilizado um refletor para detetar as posições dos tubos vazias numa calha. São dadas orientações para o funcionamento do Fluent ID por uma interface gráfica no monitor com ecrã tátil.

A radiação laser do leitor de código de barras é um raio de baixa potência, colimado no espectro visível, com as seguintes propriedades:

- Comprimento de onda: 655 nm
- Duração do impulso: 150 µs
- Potência máxima da saída de energia: 1,0 mW

4.5.1.1 Calhas de tubos compatíveis com o Fluent ID

Cada uma das calhas de tubos do Fluent ID destina-se a um tipo de tubos:

- Calha com 32 posições para tubos com 10 mm de diâmetro
- Calha com 32 posições para tubos com 13 mm de diâmetro
- Calha com 26 posições para tubos com 16 mm de diâmetro
- Calha com 32 posições para tubos Eppendorf Safe-Lock de 2 ml



É possível utilizar fichas opcionais para bloquear duas posições de uma calha de 26 posições, de forma a utilizá-la como uma calha de 24 posições, permitindo a pipetagem paralela de tubos em múltiplos de oito.





4.5.2 Manipulador FCA

Vista geral



Fig. 28: Manipulador FCA

O manipulador FCA é uma opção para o FCA configurado com adaptadores DiTi que, para além da pipetagem, permite que o FCA execute alguns movimentos de material laboratorial. O FCA consegue agarrar e soltar automaticamente os dedos do manipulador FAC durante a execução.

Dedos do manipulador FCA



Fig. 29: Dedos do manipulador FCA

Os dedos do manipulador FCA têm de ser substituídos após 2 anos ou 20000 ciclos de utilização (considera-se um ciclo o processo de agarrar, utilizar e estacionar). Os ciclos são monitorizados por um contador definido no software Fluent Control.

4 - Descrição do funcionamento Opções e dispositivos



Recetáculo da estação de encaixe do manipulador FCA



Fig. 30: Recetáculo da estação de encaixe do manipulador FCA

O recetáculo da estação de encaixe do manipulador FCA é utilizado para armazenar os dedos do manipulador FCA. Pode ser montado num recetáculo de microplaca padrão num segmento da mesa de trabalho.



4.5.3 Mistura e perfuração

A Estação de trabalho de mistura e perfuração Fluent foi concebida para aplicações que transferem líquido de e para tubos de amostras com tampas de borracha, sem que tenha de as remover, através da perfuração das respetivas tampas.



Os septos dos tubos apenas podem ser perfurados uma vez. A perfuração repetida do mesmo tubo não é suportada.

6

Os tubos BD Vacutainer® com tampa Hemogard e os tubos Greiner Vacuette® com tampa de puxar sem rebordos e tampa de rosca de segurança foram testados quanto à perfuração. A perfuração de tubos com tampões de borracha não é suportada.



As pontas de perfuração devem ser mudadas regularmente. A utilização foi verificada para 20 000 perfurações/pontas.

A Estação de trabalho de mistura e perfuração Fluent é configurada com até 2 FCA líquidas, uma estação de lavagem profunda e até 4 rotadores de tubos, dependendo do tamanho da unidade de base Fluent. Os Rotadores de tubos podem ser integrados em qualquer tamanho de unidade de base Fluent e suportam leitura de códigos de barras dos tubos, mistura das amostras, perfuração e alíquotas. Para obter mais informações sobre os tipos de tubos suportados, consulte a secção "Calhas do rotador de tubos" [▶ 78].

O curso de trabalho pode ser dividido nos seguintes passos:

- 1. Leitura de código de barras durante o carregamento do tubo
- 2. Mistura dos conteúdos dos tubos
- Perfuração e manipulação de líquidos com o FCA no Rotador de tubos com Pontas de perfuração
- 4. Lavagem e descontaminação das Pontas de perfuração na estação de lavagem profunda e cartuchos de descontaminação
- 5. Repetição do passo 2 e continuação
- 6. Recomenda-se uma lavagem prolongada da estação de lavagem, incluindo os resíduos centrais e as unidades de limpeza frontal e traseira, no final do script.

4.5.3.1 Rotador de tubos

O objetivo principal do Módulo do rotador de tubos é misturar o conteúdo líquido dos tubos e funcionar como um carregador para ações de perfuração e pipetagem. Um único Rotador de tubos tem a capacidade de 5 Calhas do rotador de tubos com 24 tubos cada (isto é, capacidade total de 120 tubos).

O dispositivo contém os seguintes subcomponentes:

 Um Leitor de código de barras de tubos integrado para ler os códigos de barras das amostras durante o carregamento



- Uma estação de lavagem profunda opcional com cartuchos profundos para descontaminação de Pontas de perfuração e um suporte de tubo de erro. O suporte de tubo de erro pode ser utilizado para guardar amostras no caso de erros de perfuração. A estação de lavagem é colocada próxima do tambor rotativo.
- Um tambor rotativo com um fixador do tubo que acomoda até cinco Calhas do rotador de tubos. O tambor realiza a mistura da amostra através de rotação ou oscilação de 360° em diferentes ângulos e velocidades. O fixador (tampa) suporta o processo de perfuração.
- O Rotador de tubos com Calhas do rotador de tubos suporta a deteção capacitiva do nível do líquido antes e após a aspiração, bem como após a distribuição de líquido (verificação da chegada do líquido) através de tubos fechados (configuração opcional).
- O Rotador de tubos é instalado pelo TASS e não deve ser movido pelo operador principal nem pelo utilizador.

4.5.3.2 Calhas do rotador de tubos

As calhas do rotador de tubos foram concebidas para uma utilização no Rotador de tubos e para suportarem a funcionalidade de perfuração. Existem diferentes tipos de calhas de tubos destinadas a acomodar os tipos de tubos suportados para perfuração:

- 13 x 75 mm Calha do rotador de tubos BD, 24 posições do tubo
- 13 x 100 mm Calha do rotador de tubos BD, 24 posições do tubo
- 13 x 75 mm Calha do rotador de tubos Greiner, 24 posições do tubo
- 13 x 100 mm Calha do rotador de tubos Greiner, 24 posições do tubo
- 16 x 100 mm Calha do rotador de tubos, 24 posições do tubo

A distribuição múltipla só é suportada para tubos Greiner Vacuette® com uma tampa de puxar sem rebordos e tampa de rosca de segurança.



Os tubos com tampões de borracha não são compatíveis e não podem ser perfurados.

Tab. 19: Compatibilidade do tubo e da calha

Linha de produtos	Tubo		Calha		
	Diâmetro [mm]	Comprim ento [mm]	Definição do material laboratorial	Calha compatível	Cor da ponte
Greiner Vacuette	13	100	13 x 100 mm Greiner Vacuette com septo	1 x 24 13 x 100 mm Calha do rotador de tubos Greiner	cinzent o
	13	75	13 x 75 mm Greiner Vacuette com septo	1 x 24 13 x 75 mm Calha do rotador de tubos Greiner	
	16	100	16 x 100 mm Greiner Vacuette com septo	1 x 24 16 x 100 mm Calha do rotador de tubos	preto

⁶



BD Vacutainer	13	100	13 x 100 mm BD Vacutainer com septo	1 x 24 13 x 100 mm Calha do rotador de tubos BD	branco
	13	75	13 x 75 mm BD Vacutainer com septo	1 x 24 13 x 75 mm Calha do rotador de tubos BD	
	16	100	16 x 100 mm BD Vacutainer com septo	1 x 24 16 x 100 mm Calha do rotador de tubos	preto

4.5.3.3 Proteção da ponta de perfuração



Fig. 31: Proteção da ponta de perfuração

A proteção da ponta de perfuração é uma tampa utilizada para cobrir o ápice afiado das pontas de perfuração durante a substituição das pontas e a resolução de problemas. Protege o utilizador contra ferimentos e as pontas contra danos.



A proteção da ponta de perfuração é de utilização única. Após a utilização, todas as proteções de pontas de perfuração têm de ser descartadas num recipiente para resíduos biológicos.





4.5.3.4 Ferramenta de remoção de pontas de perfuração

Fig. 32: Ferramenta de remoção de pontas de perfuração

A ferramenta de remoção de pontas de perfuração é utilizada para recolher uma ponta de perfuração presa num tubo que não pode ser recolhida através dos comandos do software.

4.5.4 Frida Reader



Fig. 33: Frida Reader



O Frida Reader destina-se à quantificação e normalização automatizadas dos ácidos nucleicos. As amostras a medir com o Frida Reader têm de ser arrefecidas a 4 °C para que a evaporação de amostras não comprometa os resultados das medições.

As vibrações podem causar resultados errados!

A vibração da gota da amostra pode causar resultados de medição incorretos e comprometer a segurança ou a condição clínica da amostra do paciente.

- Um pavimento estável é um pré-requisito para uma localização de instalação adequada.
- Durante as medições do Frida Reader não são permitidas fontes de vibração internas ou externas próximas.
- Evite fontes com a frequência de ressonância. Devem ser evitadas especificamente vibrações de cerca de 36 Hz (2160 rpm) e cerca de 42 Hz (2520 rpm), uma vez que estas são frequências de ressonância de uma gota suspensa.

A iluminação da divisão pode causar resultados errados!

A iluminação da divisão acima do módulo pode interferir com a medição, causar resultados de medição incorretos e comprometer a segurança ou a condição clínica da amostra do paciente.

• O sistema robótico tem de ter uma tampa não transparente e um painel frontal e traseiro para evitar a luz ambiente na posição de medição do Frida Reader.

4.5.5 Separador de fases

O separador de fases foi concebido para detetar fases de separação entre líquidos de viscosidade diferente. Como tal, é independente de a fase de separação ser visível a partir do exterior do material laboratorial. O separador de fases pode ser utilizado em aplicações que requerem a transferência limpa de uma fase líquida de um material laboratorial de origem para um de destino.

O que se segue representa um fluxo de trabalho típico:

- Centrifugação do material laboratorial de origem com uma mistura de líquidos para gerar uma fase distinta entre líquidos.
- Os líquidos têm de diferir na sua viscosidade para permitir a formação de fases durante a centrifugação.
- Carregue os tubos sem tampa para a plataforma do Fluent e utilize um leitor de códigos de barras (por exemplo, Fluent ID) para uma rastreabilidade total. Proceder com cuidado para não perturbar a camada entre as fases/frações líquidas durante o carregamento.
- Iniciar o protocolo definido para a separação das frações. A função de separador de fases do Air FCA deteta a fase entre os líquidos e inicia a transferência da fração de interesse para o material laboratorial de destino.



É possível extrair mais do que uma fase do material laboratorial de origem. A remoção de, pelo menos, parte da fase superior é necessária para evitar derrames de líquido (extravasamento de material laboratorial) na plataforma e a potencial contaminação do canal de pipetagem acima da ponta descartável durante a deteção de fases.

Para obter as especificações técnicas, consultar o manual de referência. Para mais informações sobre o software FluentControl, consulte o manual do software de aplicação. Consulte o "Documentos de referência" [> 9].

As fases líquidas têm de estar claramente separadas. Para a separação do sangue total do plasma e células sanguíneas, a qualidade e o pré-tratamento das amostras são cruciais. Os parâmetros que podem ter um efeito na deteção de fases em amostras de sangue são a qualidade das amostras (lipemia, hemólise), o tempo de conservação, a temperatura de conservação, as condições de

transporte, as condições de centrifugação (tempo, rcf, temperatura, rampa, tipo de rotor), a distorção da fase após a centrifugação, etc.

Para obter os melhores resultados de separação de fases, as amostras de sangue devem ser processadas o mais rapidamente possível após a colheita. As condições de tratamento e conservação das amostras devem seguir as recomendações do fabricante do tubo específico.

A centrifugação a 2500 rcf durante 10 minutos à temperatura ambiente com uma rampa de desaceleração resulta numa separação de fase limpa para amostras de plasma (condições de teste internas Tecan).

Para evitar o derrame de líquido dos tubos durante a aspiração e a deteção de fases, os tubos não devem ser enchidos até à extremidade. Por norma, a deteção de fases requer um movimento descendente rápido no tubo, combinado com uma velocidade de aspiração lenta, o que leva a um aumento do nível de líquido durante a deteção.

Para tubos com um volume de enchimento elevado, recomenda-se a remoção de líquido da parte superior antes de se iniciar a deteção de fases.



4.5.6 Resolvex i300



Fig. 34: Módulo da mesa de trabalho Resolvex i300

O Resolvex i300 é um módulo que pode ser integrado na plataforma Fluent. A plataforma Fluent pode, então, manusear líquidos e transferir material laboratorial a partir de uma mesa de trabalho diretamente para o módulo Resolvex i300, dentro da mesma estrutura de software.

O Resolvex i300 é um instrumento de laboratório que permite automatizar vários processos laboratoriais utilizando as seguintes funções principais:

- Pressurização das colunas do filtro (total ou parcialmente carregadas) com ar ou nitrogénio
- Empilhamento e desempilhamento de placas de filtro em placas de colheita com um elevador de material laboratorial integrado
- Proteção das amostras contra contaminação cruzada durante o empilhamento e desempilhamento de material laboratorial de filtração e colheita utilizando um protetor contra pingas integrado
- Dispensação de líquidos para material laboratorial de filtração e colheita (opcional)
- Evaporação de líquidos do material laboratorial de colheita com gás aquecido, como ar ou nitrogénio (opcional)



Os fluxos de trabalho de aplicação típicos do Resolvex i300 incluem os seguintes processos laboratoriais:

- Processos de extração em fase sólida (SPE) para fluxos de trabalho de preparação de amostras de espectrometria de massa
- Extração de líquidos suportada (SLE)
- Outros fluxos de trabalho de filtração, purificação ou concentração que utilizam pressão positiva

Para obter mais informações e antes de executar qualquer trabalho com o Resolvex i300, consulte o Manual de operação do Resolvex i300. Consulte o "Documentos de referência" [▶ 9].



5 - Elementos de controlo Elementos de comando

5 Elementos de controlo

5.1 Elementos de comando



As alavancas de grampo bloqueiam e desbloqueiam os segmentos.

O ecrã tátil apresenta os métodos e as descrições, permitindo ao operador controlar o instrumento.



5.2 Interface do utilizador



Fig. 36: Interface do utilizador do FluentControl

- A Caminho de navegação B Área de trabalho
- C Botões do mostrador/de opção/de ação

Através da interface do utilizador do FluentControl, o operador tem acesso às execuções do método para a operação e os cuidados do sistema.

5.2.1 Caminho de navegação

Utilize o caminho de navegação para compreender e percorrer a hierarquia do FluentControl.

Botão	Nome	Função
$ \Delta\rangle$	Home	Prima para regressar à página inicial.
System Care 🛞	Navigation pane	Apresenta as seleções atuais e as anteriores.
<	Menu expander	Prima o botão do expansor do menu para revelar opções como os controlos das luzes e para mudar de operador.

Tab. 20: Botões do caminho de navegação



5.2.2 Área de trabalho

Aceda a métodos e descrições através da área de trabalho da interface do utilizador. Aqui são também apresentados detalhes acerca do estado da execução do método.

Tab. 21: Botões da área de trabalho

Botão	Nome	Função
	Run	Prima para iniciar o método selecionado.
+	Add	Prima para adicionar mais métodos à sua lista de início rápido.
Assay 1 urknown	Selected method	O método atualmente selecionado que será executado quando pressionar o botão Continue.
Assay 3 unknown	Available method	Um método que pode ser selecionado se clicar no mesmo.
Start of Day Due: is not executed	Botão Quick start	Prima para iniciar imediatamente o método selecionado.

Tab. 22: Mostrador da área de trabalho

Mostrador	Função do mostrador
00:16:52 10:09 10:43 Star Time Tai Time	Apresenta o estado e o tempo restante da execução do método.
Assay 1 is ready to be started.	Descrição do método atualmente selecionado ou informações adicionais sobre a ação atual.



5.2.3 Botões do mostrador, de opção e de ação

Tab. 23: Botões do mostrador, de opção e de ação

Botão	Nome	Função
√ 0k	Ok	Prima para confirmar.
× Cancel	Cancel	Prima para cancelar.
Continue	Continue	Prima para continuar.
II Pause	Pause	Prima para pedir uma pausa na execução no final da ação atual.
Stop	Stop	Prima para parar uma execução imediatamente, mesmo a meio da ação atual. Se possível, o sistema irá oferecer a possibilidade de restaurar ou de continuar a execução.
THE Remove	Remove	Prima para remover o método da vista de início rápido.
View Mode	View mode	Prima para alternar entre a vista da lista e as vistas de início rápido.
Sort by	Sort by	Prima para alternar a apresentação da execução do método entre a ordenação alfabética e pela mais recente.

5.2.4 Botões de recuperação do método

Tab. 24: Botões do mostrador, de opção e de ação

Botão	Nome	Função
Discard	Discard	Prima para eliminar o estado de um método recuperado.



Botão	Nome	Função
\rightarrow	_	Prima para passar para o ecrã seguinte.
Recovery Point	Recovery Point	Prima para regressar ao ecrã anterior ("Recovery Point").
Run Recovery	Run Recovery	Prima para continuar a execução.

5.2.5 Botões DeckCheck

Tab. 25: Botões DeckCheck

Botão	Nome	Função
Left	Câmara esquerda	Apresenta a imagem capturada pela câmara esquerda (apenas no Fluent 780/1080). Um ponto de exclamação surge no ícone se tiver sido observada uma discrepância de disposição com esta câmara.
Center	Câmara central	Apresenta a imagem capturada pela câmara central de vista geral. Um ponto de exclamação surge no ícone se tiver sido observada uma discrepância de disposição com esta câmara.
Right	Câmara direita	Apresenta a imagem capturada pela câmara direita (apenas no Fluent 780/1080). Um ponto de exclamação surge no ícone se tiver sido observada uma discrepância de disposição com esta câmara.
Pause Alternate	Pausar alternância	O ecrã alterna entre imagens ao vivo e de referência: Prima este botão quando está a ser apresentada a imagem ao vivo ou de referência para manter essa imagem estática.



Botão	Nome	Função
Resume Alternate	Retomar alternância	A imagem está estática: Prima para retomar a alternância entre as imagens ao vivo e de referência.
Check	Verificar	Ativa uma nova verificação do sistema – por exemplo quando foram feitas algumas correções. Será solicitado que feche a porta. Para um sistema de 3 braços, o braço central deve mover-se: Se a porta não estiver fechada, a verificação será realizada, mas o braço central irá bloquear uma câmara.
Ignore & Continue	Ignorar e continuar	Apenas é apresentado se estiver configurado para o respetivo comando no método. Permite que sejam ignoradas quaisquer discrepâncias realçadas, e a execução do guia irá continuar.
Continue	Continuar	É apresentado quando todas as discrepâncias tiverem sido resolvidas ou se o sistema não tiver encontrado nenhuma discrepância e a opção Show always tiver sido selecionada para o comando. Tal pode permitir que alterações de cor subtis que o sistema não reconheceu sejam visíveis a olho nu.

Tab. 26: Visores

Visor	Descrição	Função
	Imagem de referência	A imagem de referência é armazenada no comando da guia com a disposição da mesa de trabalho pretendida.
LIVE	Imagem ao vivo	A imagem ao vivo capturada pelas câmaras durante a execução da guia.
	Discrepância (diferença em relação à imagem de referência)	Os quadrados vermelhos assinalam as áreas onde foram encontradas discrepâncias entre as imagens ao vivo e de referência. A área assinalada pode incluir mais do que um erro.





5.3 Sinais de erro e estado do instrumento

Fig. 37: Lâmpadas de estado

- A Lâmpada de estado da energia B
- Lâmpada de estado superior



As lâmpadas de estado indicam o estado do instrumento através de luzes de diferentes cores, fixas ou intermitentes. A lâmpada de estado superior está ativa apenas quando o software está a ser executado.

Tab. 27: Sinais luminosos das lâmpadas de estado

Sinal	Cor	Modo	Estado do instrumento
	_	desligado	O instrumento encontra-se desligado (desconectado da fonte de alimentação).
	branco	"pulsação"	O instrumento encontra-se ligado (software de controlo conectado, módulos ainda não iniciados).
	branco (lâmpada de energia apenas)	contínuo	Instrumento no estado "ligado" (software de controlo ainda não conectado).



Sinal	Cor	Modo	Estado do instrumento
	esquema de	"pulsação"	Modo de inatividade
••••	cores da interface do utilizador do FluentControl		Todos os módulos estão inicializados; o instrumento está pronto para executar um método.
			Após cerca de uma hora no modo de inatividade, o instrumento irá passar para o modo de standby.
			Modo de standby
			Todos os eixos estão travados. Os braços não se encontram no ZeroG e não podem ser movidos manualmente. Para ativar o instrumento, execute um método ou solicite ao operador principal para selecionar a ferramenta de deslocamento para o modo ZeroG.
	amarelo	contínuo	Modo de configuração
			O instrumento "aprende" as posições.
			Neste modo, o utilizador pode deslocar manualmente os braços robóticos.
	verde	contínuo	Está em execução um método (guia ou processo).
			Este é o modo de "produção" regular.
	vermelho	intermitente	Estado de erro
			O ecrã do computador de controlo ou o ecrã tátil apresenta uma mensagem de erro.
	a cor é	intermitente	Solicitação do utilizador
	configurável pelo utilizador		O sistema aguarda uma interação do utilizador.
	verde	intermitente	Paragem ativa
			Esta é uma pausa intencional desencadeada pelo controlador de tempo de execução ou pela abertura de um painel de segurança.
			O instrumento entra em pausa para permitir a interação do utilizador com a mesa de trabalho. O operador pode retomar o método.



5.4 LED de estado do Fluent ID



Fig. 38: LED do Fluent ID

Os LED do Fluent ID assinalam os seguintes estados:

Tab. 28: LED do Fluent ID

Sinal	Cor	Modo	Estado do instrumento
	_	desligado	O Fluent ID está inativo.
	branco	contínuo	O Fluent ID está ligado (mas ainda não iniciado).
	azul ou cor personalizada	intermitente	Pronto para carregar ou descarregar a calha.
	verde	contínuo	Códigos de barras lidos com sucesso. Calha supervisionada. Não descarregue, uma vez que essa ação irá interromper a execução.
	vermelho	intermitente	Estado de erro A mensagem de erro e a ação necessária são apresentadas no ecrã tátil.



6 Funcionamento

6.1 Instruções de segurança deste capítulo

Resultado errado ou contaminação do instrumento!

Poderão ocorrer resultados errados ou a contaminação do instrumento se a qualificação da instalação e a qualificação operacional não tiverem sido realizadas ou se os procedimentos operativos indicados neste manual não forem seguidos.

- Os registos da qualificação da instalação e da qualificação operacional estão disponíveis e são conhecidos.
- Os métodos e os processos, incluindo os parâmetros de pipetagem, devem ser validados pelo operador principal.
- A deteção do nível de líquido, em conjunto com as aplicações de perfuração para o FCA e o Air FCA devem ser validados pelo operador principal.
- Para aplicações de sangue total com pontas de perfuração, rotador de tubos e soluções concentradas de descontaminação corrosivas, recomenda-se a substituição de ambos os conectores da estação de lavagem (na estação de lavagem e no recipiente de líquido do sistema) a cada três meses para evitar obstruções e desgaste dos conectores. Os líquidos potencialmente corrosivos têm de ser neutralizados antes da drenagem, ou a estação de lavagem tem de ser purgada com líquidos neutros, como água.
- O operador deve receber formação relativamente aos procedimentos operativos, métodos e processos.



Contaminação biológica e química do utilizador!

Os dedos do manipulador FCA danificados podem perder placas. As placas que caíram podem provocar contaminação por substâncias perigosas.

• Verifique os dedos do manipulador FCA após um embate.

▲ CUIDADO

Extremidades e pontas afiadas!

As Pontas de perfuração da Estação de trabalho de mistura e perfuração Fluent possuem cabeças pontiagudas e extremidades afiadas que podem provocar ferimentos.

- Ao carregar o instrumento, desloque o FCA para uma posição segura através de um comando de software.
- Quando ocorrer um erro, cubra as pontas de perfuração com proteções de pontas de perfuração e desloque o FCA manualmente para uma posição segura. Consulte a secção "Proteção da ponta de perfuração" [▶ 79].

▲ CUIDADO

Contaminação biológica do sistema!

Na Estação de trabalho de mistura e perfuração Fluent, o sangue pode contaminar as tampas dos tubos.

- Manuseie os tubos com cuidado.
- Use equipamento de proteção.

AVISO

Avaria devido a líquidos corrosivos!

A drenagem de líquidos corrosivos concentrados, como lixívia a 2%, através da estação de lavagem e dos tubos pode provocar avaria dos conectores da estação de lavagem.

 Os líquidos potencialmente corrosivos têm de ser neutralizados antes da drenagem, ou a estação de lavagem tem de ser purgada com líquidos neutros, como água.

6.2 Modos de funcionamento

O Fluent pode trabalhar em três modos de funcionamento diferentes:

Operador

Modo de funcionamento de rotina

- Modo de funcionamento normal, em que a aplicação ou as tarefas de cuidados de rotina do sistema são executadas.
- O Fluent é monitorizado pelo controlador runtime do software do FluentControl.

FSE



Operador
principalModo de definição do método• Este modo de funcionamento é utilizado para realizar tarefas especiais, tais

como o ajuste para configurar o método.

Modo de assistência

- Este modo de funcionamento é utilizado para realizar tarefas especiais, tais como os testes para assegurar a operacionalidade.
- É necessário um certificado de assistência para executar este modo.

O braço embate com objetos na mesa de trabalho

Os braços do Fluent podem ser movidos manualmente. Certifique-se de que os movimentos manuais dos braços são efetuados com suavidade, mantendo o braço fixo. Não deixe os braços embaterem com objetos sólidos, incluindo o batente mecânico do alcance do braço

6.3 Colocação em funcionamento

6.3.1 Ligar o instrumento

Para ligar o instrumento, proceda do seguinte modo:

1. Ligue a energia no interruptor de energia (A) na traseira da fonte de alimentação externa.





Quando o instrumento ligar, a lâmpada de energia irá acender a azul. Consulte a secção "Sinais de erro e estado do instrumento" [▶ 91].

Se a lâmpada de estado não acender, ligue o PC ou contacte o operador principal.

 Inicie o software do FluentControl. Consulte a secção "Iniciar o FluentControl" [> 97].

6.3.2 Iniciar o FluentControl

- ✓ Os procedimentos operativos devem estar disponíveis e ser conhecidos.
- Os registos da qualificação da instalação e da qualificação operacional estão disponíveis e são conhecidos.
- ✓ Os cuidados do sistema foram realizados.
- ✓ O instrumento está ligado.
- Inicie o software com Start > All programs > Tecan > FluentControl. Após alguns segundos, o ecrã Start é apresentado.

6.3.3 Início de sessão do utilizador

Para iniciar sessão no FluentControl, proceda do seguinte modo:

- ✓ O software Fluent Dx está instalado.
- ✓ O FluentControl foi iniciado.
- ✓ A Gestão do utilizador está ativada no FluentControl e o processo foi definido.
- 1. Selecione o perfil do utilizador atribuído (A).



2. Introduza a palavra-passe no teclado (B).



3. Prima OK (C).



Após o início de sessão, o instrumento é automaticamente inicializado.

6.3.4 Colocar segmentos

Para colocar segmentos, proceder do seguinte modo:

- ✓ Todos os segmentos, transportadores, opções e dispositivos têm de ser colocados de acordo com o método selecionado.
- ✓ Os segmentos estão limpos e em perfeitas condições.
- ✓ Os segmentos estão colocados na posição da grelha correspondente.
- 1. Baixar o segmento para a parte traseira da plataforma.
- 2. Alinhar a extremidade traseira com a tampa do canal traseiro ou a extensão do instrumento.





3. Baixar cuidadosamente a parte dianteira do segmento (A).

4. Rodar a alavanca de grampo da esquerda para a direita, para a posição fechada. Consulte a secção "Verificar o segmento" [▶ 206].

6.3.5 Remover segmentos



Os segmentos do Fluent ID não foram concebidos para serem removidos! Estes estão conectados diretamente à parte eletrónica do instrumento. A conexão à parte eletrónica pode ser realizada apenas por um FSE qualificado. Consulte a secção Apoio ao cliente.

Para remover segmentos, proceda do seguinte modo:

- ✓ Todos os reagentes, amostras, racks, calhas e placas foram removidos do segmento.
- ✓ Não se encontra nada colocado no segmento.
- 1. Rode a alavanca de grampo (A) da direita para a esquerda, para a posição aberta.





O segmento encontra-se desbloqueado e a marca amarela na alavanca de grampo está visível.



- 2. Empurre o segmento para a frente, aproximadamente 4 mm.
- 3. Levante o segmento (B) na parte da frente.



4. Armazene o segmento num local limpo e seco, para evitar danos.

6.3.6 Carregar as calhas standard

AVISO

Danos devido ao carregamento ou descarregamento inadequado

Danos nas calhas e nos pinos.

- Alinhe a calha horizontalmente com a plataforma.
- Sustenha a parte da frente da calha com uma mão.
- Ao descarregar, certifique-se de que a calha liberta todos os pinos antes de a levantar.

Para carregar as calhas, proceda do seguinte modo:





1. Alinhe a calha com a respetiva posição da grelha (A).

- 2. Empurre a calha para a posição de paragem.
- Certifique-se de que a calha bloqueia o segmento de forma segura. Pode senti-lo nos últimos milímetros antes de a calha tocar na posição de paragem.



Para descarregar as calhas, proceda do seguinte modo:

- 1. Puxe a calha horizontalmente ao nível da mesa de trabalho até esta ser totalmente removida da área de carregamento.
- 2. Sustenha a parte da frente da calha com uma mão.
- 3. Certifique-se de que a calha liberta todos os pinos antes de a levantar.



6.3.7 Verificar a disposição da mesa de trabalho

Certifique-se de que os carregadores, o material laboratorial e os dispositivos instalados na mesa de trabalho correspondem à disposição da mesa de trabalho definida para o método.

AVISO

Danos no equipamento!

O posicionamento incorreto do segmento e do material laboratorial na mesa de trabalho pode causar o embate dos braços.

- Certifique-se sempre de que a configuração da plataforma física e o material laboratorial carregado correspondem à configuração da mesa de trabalho do FluentControl.
- Certifique-se sempre de que o material laboratorial é instalado corretamente nos recetáculos. Consulte "Posição do material laboratorial" [> 180].

AVISO

O campo magnético cria interferência!

Um campo magnético forte (polo norte para cima) na posição de aspiração pode interferir com o sensor da presença de pontas e pode levar a erros imprevistos (por ex., **perda de DiTi**).

• Certifique-se de que não se encontra nenhum íman forte colocado numa posição SBS adjacente à posição de aspiração.



Os segmentos do Fluent ID só podem ser removidos pelo TASS, devido à conexão ao quadro elétrico por baixo da mesa de trabalho.

- ✓ O método deve ser preparado pelo operador principal.
- ✓ Os consumíveis são consistentes com os consumíveis definidos no método.
- As calhas de tubos do Fluent ID devem ser carregadas apenas após o início do método, quando solicitado no ecrã tátil.
- 1. Siga as instruções apresentadas no ecrã tátil.



A ilustração apresenta um exemplo de uma instrução (A) apresentada no ecrã tátil:



6.4 Antes de iniciar um método

A lista de verificação seguinte tem de ser concluída antes de iniciar um método.

Instrumento/ Componente	Tarefa	Referência/Atividades
Validação do processo	Antes de iniciar uma execução de produção, certifique-se de que o método que está a selecionar foi validado.	Contacte o operador principal para obter mais informações.
Ecrã táctil	Siga as instruções do ecrã tátil. AVISO! As instruções disponibilizadas pelo operador principal têm de ser cumpridas rigorosamente.	
	Caso não sejam apresentadas quaisquer instruções, siga a lista de tarefas seguinte.	



Instrumento/ Componente	Tarefa	Referência/Atividades
Segmentos, transportadores, opções e dispositivos	Certifique-se de que todos os segmentos, transportadores, opções e dispositivos estão instalados e seguros. Certifique-se de que se encontram na plataforma apenas os objetos destinados a utilizar no método. Certifique-se de que a execução de teste é concluída com sucesso.	Se a execução de teste falhar, contacte o operador principal para realizar a execução de teste novamente.
Amostras e reagentes	Certifique-se de que todas as amostras, reagentes e material laboratorial são carregados corretamente. AVISO! A leitura de códigos de barras acontece apenas após o início do método. Certifique-se de que a plataforma do Fluent ID não tem calhas antes do início do método. As calhas só podem ser carregadas quando o pedido é apresentado no ecrã tátil.	
Tubagem de desperdícios (apenas nos sistemas de líquido)	Certifique-se de que a tubagem de desperdícios está posicionada corretamente.	Inspecione visualmente a tubagem de desperdícios para se certificar de que não se encontra dobrada ou esmagada.
Sistema de lavagem (apenas nos sistemas de líquido)	Certifique-se de que o líquido do sistema e o recipiente de desperdícios se encontram corretamente ligados.	Consulte a secção "Verificar a tubagem do recipiente do líquido do sistema e do recipiente para resíduos" [▶ 106].





Instrumento/ Componente	Tarefa	Referência/Atividades
Sistema de lavagem (apenas nos sistemas de líquido)	Certifique-se de que o recipiente do líquido do sistema está cheio até ao nível correto. Certifique-se de que o recipiente de desperdícios está vazio.	Consulte a secção "Conectar o recipiente do líquido do sistema e o recipiente para resíduos" [▶ 158].
Sistema de lavagem (apenas nos sistemas de líquido)	Certifique-se de que é utilizado o líquido do sistema correto, conforme definido no método.	_
Desperdícios de pontas descartáveis e unidade da estação de lavagem	Certifique-se de que os desperdícios de pontas descartáveis e a unidade da estação de lavagem se encontram limpos.	Consulte a secção "Limpar os resíduos de pontas descartáveis e a unidade da estação de lavagem" [▶ 152].
	Certifique-se de que as tampas das calhas de resíduos para a contenção de aerossóis estão montadas.	_
Pontas descartáveis	Certifique-se de estão carregadas as pontas descartáveis corretas. Certifique-se de que a unidade de desperdícios de pontas está vazia.	
Pontas fixas	Certifique-se de que as pontas fixas estão limpas e sem danos.	Inspecione visualmente as pontas fixas para se certificar de que se encontram limpas. Inspecione visualmente as pontas fixas com um espelho de dentista, para se certificar de que o revestimento está intacto.
Plataforma	Certifique-se de que os transportadores, o material laboratorial e os dispositivos instalados na plataforma correspondem à disposição da plataforma definida para o método.	Consulte a secção "Verificar a disposição da mesa de trabalho" [▶ 102].



Instrumento/ Componente	Tarefa	Referência/Atividades
Material laboratorial	Certifique-se de que todo o material laboratorial está posicionado de forma segura. Se as microplacas apresentarem um movimento lateral, certifique-se de que os posicionadores de material laboratorial estão corretos.	Consulte a secção "Posição do material laboratorial" [▶ 180].
Rotador de tubos	Certifique-se de que não há pinos de posicionamento ou pinos de bloqueio em falta no Rotador de tubos.	Consulte a secção "Substituir os pinos de bloqueio e os pinos de posicionamento" [▶ 208]

6.4.1 Verificar a tubagem do recipiente do líquido do sistema e do recipiente para resíduos

Contaminação da amostra!

No caso de uma configuração dupla do Liquid FCA, podem ser utilizados diferentes tipos de líquidos do sistema para cada braço. Ligar o recipiente do líquido do sistema errado a um braço pode causar a contaminação da amostra.

- Identifique cada recipiente do líquido do sistema com o nome do líquido do sistema correspondente.
- ✓ O sistema de lavagem deve ser instalado corretamente.
- 1. Verifique se o tubo (B) está corretamente ligado ao recipiente do líquido do sistema (D).





2. Caso esteja presente, verifique se o sensor de deteção de líquido (C) está corretamente ligado ao recipiente do líquido do sistema (D).

6.4.2 Verificar a tubagem do recipiente para resíduos

1. Verifique se o tubo (A) está ligado ao recipiente para resíduos (D).



- 2. Verifique se os tubos (C) e (D) se encontram corretamente ligados ao recipiente para resíduos.
- 3. Caso esteja presente, verifique se o sensor de deteção de líquido (E) está corretamente ligado ao recipiente para resíduos.



4. Enrosque a tampa (F).



6.5 Executar um método

Um método é uma recolha de guias ou processos definidos no software do FluentControl. O método pode ser executado numa execução.

O operador principal escreve um método que pode ser executado como se segue.

AVISO

Danos no instrumento!

Poderão ocorrer danos no instrumento se a mesa de trabalho não estiver corretamente configurada ou se o software for utilizado ou operado incorretamente.

- Certifique-se de que todos os dispositivos de segurança estão instalados e funcionais.
- Certifique-se de que os carregadores, o material laboratorial e os dispositivos instalados na mesa de trabalho correspondem à disposição da mesa de trabalho definida para o método.
- Certifique-se de que se encontram na mesa de trabalho apenas os objetos destinados a utilizar no método.

6.5.1 Iniciar um método

- ✓ O FluentControl foi iniciado.
- ✓ A secção Início de sessão do utilizador foi executada.
- ✓ A secção foi executada.
- 1. Selecione Method Starter (Iniciador de método) (A).






- Selecione o método (B) a executar.
 O método selecionado é realçado.
- 3. Prima OK.





4. Prima Run (Executar) (C).



- 5. Siga as instruções do ecrã tátil.
- 6. Se a sua guia incluir o DeckCheck, preste atenção a quaisquer diferenças na disposição da mesa de trabalho real em comparação com a disposição da mesa de trabalho esperada da mesa de trabalho de referência. Consulte Operação DeckCheck.
- 7. Aguarde que a execução do método termine.

O ecrã apresenta o tempo aproximado (D) até ao fim da execução do método.





- Construction
 Running

 Run finished with error(s)!

 11:59
 12:19

 Star Time
 12:19

 Star Time
 12:19

 Image: Comment Process
 Log

 More
 Process
- 8. Se for apresentada a mensagem **Run finished with error(s)!** (Execução terminada com erro[s]!), prima **Log** (Registo) para rever os erros e avisos.

9. Prima Next (Seguinte) para regressar ao ecrã Home (Início).



10. Prima Exit (Sair).





O ecrã abaixo (E) é apresentado quando a execução do método termina.

6.5.2 Carregar e descarregar as calhas do Fluent ID



O Fluent é um produto laser de classe 1, de acordo com a norma IEC 60825-1:2014, que emite radiação laser.

O raio laser pode causar encadeamento, cegueira por ofuscamento e imagens persistentes.

Não olhe para o raio laser nem para os reflexos deste.

6.5.2.1 Carregar as calhas do Fluent ID

AVISO

Danos devido ao carregamento ou descarregamento inadequado

Danos nas calhas e nos pinos.

- Alinhe a calha horizontalmente com a plataforma.
- Sustenha a parte da frente da calha com uma mão.
- Ao descarregar, certifique-se de que a calha liberta todos os pinos antes de a levantar.
- ✓ O Fluent está equipado com um leitor de códigos de barras de tubos Fluent ID.
- ✓ Os tubos estão carregados nas calhas com as etiquetas dos códigos de barras voltadas para a esquerda.
- ✓ Todos os tubos numa calha têm o mesmo tamanho e formato. Para obter informações sobre os tipos de calhas de tubos, consulte a secção "Calhas de tubos compatíveis com o Fluent ID" [▶ 74].
- 1. Selecione e inicie o método, utilizando o ecrã tátil.

Os LED começam a piscar e a mensagem **Please load tubes** é apresentada no ecrã tátil.



Se utilizar diferentes tipos de tubos, certifique-se de que é utilizado o tipo de calha correto para cada grelha específica.

- 2. Sustenha a parte da frente da calha com uma mão.
- 3. Segure a calha horizontalmente ao nível da plataforma.
- 4. Empurre a calha para a posição de paragem.



5. Faça deslizar as calhas de tubos, uma após a outra, nas grelhas dedicadas na área de carregamento do Fluent ID.

$\left egin{array}{c} \Delta \end{array} ight angle ight angle$ Running $ ight angle$ Load tub	e runners		
		000000000000000000000000000000000000000	
	1 2 3 4	5 6	
Please load tube runner	s and press continue when	finished	Errors: 🔲 Missing Runner
			V V
			Stop Continue

Verifique se todos os códigos de barras foram lidos com sucesso.
 Os LED ficam verdes quando as calhas de tubos se encontram na posição carregada e todas as etiquetas de códigos de barras foram lidas com sucesso.





Para obter informações sobre a descrição do estado do LED do Fluent ID, consulte a secção "LED de estado do Fluent ID" [> 93].

- 7. Em caso de erro de leitura de um código de barras, descarregue a calha, corrija o problema e carregue a calha novamente.
- 8. Puxe a calha horizontalmente ao longo da plataforma, até esta ser totalmente removida.



O Fluent ID lê todos os códigos várias vezes à medida que passam pelo leitor. Para tubos pequenos e estreitos (isto é, diâmetro \leq 10 mm), reduza a velocidade do carregamento manual para permitir todas as leituras e reduzir os relatórios de erros.

00000000000000000		000000000000000000000000000000000000000	0000000000000	000000000000000000000000000000000000000	000000000000000000000000000000000000000			
1	2	3	4	5	6			

Errors: 🔲 Missing Runner 🗋 Wrong Runner ● Unreadable 单 Duplicates 오 Missing

Fig. 39: Confirmação da leitura do código de barras apresentada no ecrã tátil

Tab. 30: Significado do GUI (calha)

Quadrado (calha)	Significado	
Verde	Todos os códigos de barras dos tubos da calha lidos com sucesso.	



Quadrado (calha)	Significado
Branco com contorno vermelho	Tipo de calha errada para esta posição da grelha.
Cinzento com contorno vermelho	Calha em falta. Deve ser carregada uma calha nesta posição da grelha.

Tab. 31: Significado do GUI (posição do tubo)

Círculo (posição do tubo)	Significado
Verde	Códigos de barras lidos com sucesso.
Vermelho	Código de barras ilegível
Laranja	Código de barras duplicado
Branco com contorno vermelho	Tubo em falta. Deve ser carregado um tubo nesta posição.



Quando é utilizada a calha de tubos Safe-Lock de 2 ml, não é possível diferenciar entre os tubos em falta e os códigos de barras ilegíveis. Os tubos em falta são reportados como códigos de barras ilegíveis.

6.5.2.2 Descarregar as calhas do Fluent ID

- ✓ A execução terminou ou está em curso uma execução e os LED piscam com a mensagem Please unload tubes apresentada no ecrã tátil.
- 1. Puxe a calha horizontalmente ao longo da mesa de trabalho, até esta ser totalmente removida.

6.5.3 Carregar e descarregar as Calhas do rotador de tubos

6.5.3.1 Carregar as Calhas do rotador de tubos

Biocontaminação do sistema e/ou do utilizador!

Os tubos de amostras danificados podem implodir, dando origem ao derramamento de amostras no Rotador de tubos.

- Certifique-se de que nenhum tubo danificado é carregado no Rotador de tubos.
- ✓ O Fluent está equipado com um Rotador de tubos.
- Os tubos são carregados nas calhas do Rotador de tubos com a etiqueta de código de barras voltada para a esquerda.
- ✓ Todos os tubos numa calha têm o mesmo tamanho e formato. Para obter informações sobre os tipos de calhas de tubos, consulte a secção "Calhas do rotador de tubos" [▶ 78].
- 1. Selecione e inicie o método, utilizando o ecrã tátil.



Os LED começam a piscar e a mensagem **Please load tubes** é apresentada no ecrã tátil.

Quando utilizar diferentes tipos de tubos, certifique-se de que escolhe a calha correta de acordo com o tipo de tubo (BD ou Greiner). Certifique-se igualmente de que carrega os tubos de alturas diferentes nas calhas correspondentes: os tubos são sempre mantidos em posição pela ponte da calha à altura das respetivas tampas. As partes inferiores dos tubos têm de assentar sempre e de forma correta nas inserções dos tubos das calhas.

2. Abra a alavanca de bloqueio da calha.



3. Sustenha a parte da frente da calha com uma mão.



- 4. Segure a calha horizontalmente ao nível da plataforma.
- 5. Empurre a calha para a posição de paragem.



6. Faça deslizar as calhas do Rotador de tubos, uma após a outra, nas grelhas dedicadas no Rotador de tubos.



7. Verifique se todos os códigos de barras foram lidos com sucesso.

Os LED ficam verdes quando as calhas dos tubos são colocadas em posição e todas as etiquetas de código de barras foram lidas com sucesso. Para obter a descrição do estado do LED do Rotador de tubos, consulte a secção "LED de estado do Fluent ID" [> 93].

- 8. Em caso de erro de leitura de um código de barras, descarregue a calha, corrija o problema e carregue a calha novamente.
- 9. Feche a alavanca de bloqueio da calha.





00000000000		000000000000000000000000000000000000000	0000000000000	000000000000000000000000000000000000000	000000000000000000000000000000000000000	
1	2	3	4	5	6	
Errors	: 🗆	Missi	ng Ru	Inner		Wrong Runner Unreadable Duplicates Missing

Fig. 40: Confirmação da leitura do código de barras apresentada no ecrã tátil

Tab. 32: Significado do GUI (calha)

Quadrado (calha)	Significado
Verde	Todos os códigos de barras dos tubos da calha lidos com sucesso.
Branco com contorno vermelho	Tipo de calha errada para esta posição da grelha.
Cinzento com contorno vermelho	Calha em falta. Deve ser carregada uma calha nesta posição da grelha.

Tab. 33: Significado do GUI (posição do tubo)

Círculo (posição do tubo)	Significado
Verde	Códigos de barras lidos com sucesso.
Vermelho	Código de barras ilegível
Laranja	Código de barras duplicado
Branco com contorno vermelho	Tubo em falta. Deve ser carregado um tubo nesta posição.

6.5.3.2 Descarregar as Calhas do rotador de tubos

Não armazene Calhas do rotador de tubos carregadas com tubos que não cumprem as condições de operação de manipulação de líquidos. Consulte a secção "Condições ambientais" [> 45].

- A execução terminou ou está em curso uma execução e os LED piscam com a mensagem Please unload tubes apresentada no ecrã tátil.
- ✓ O Rotador de tubos está na posição inicial horizontal.



1. Abra a alavanca de bloqueio da calha.



2. Puxe a calha horizontalmente ao longo da mesa de trabalho, até esta ser totalmente removida.



Os segmentos do Rotador de tubos apenas podem ser removidos pelo TASS, devido à conexão ao quadro elétrico por baixo da mesa de trabalho.

- O método deve ser preparado pelo operador principal.
- Os consumíveis são consistentes com os consumíveis definidos no método.
- As Calhas do rotador de tubos devem ser carregadas apenas após o início do método, quando solicitado no ecrã tátil.

6.5.4 Repor os erros

Se for apresentada uma mensagem, proceda do seguinte modo:

Mensagem

- Verifique a função do mostrador, a função do botão ou a mensagem de erro. Consulte as secções "Área de trabalho" [▶ 87] e "Botões de recuperação do método" [▶ 88].
- 2. Siga as instruções deste manual e do ecrã tátil para corrigir o erro.
- 3. Continue a execução do método. Consulte a secção "Botões do mostrador, de opção e de ação" [▶ 88].

Se a lâmpada de estado acender ou mudar de cor, proceda do seguinte modo:

Lâmpada de estado

- Verifique o estado do instrumento. Consulte a secção "Sinais de erro e estado do instrumento" [▶ 91].
- Se o Fluent estiver equipado com um leitor de códigos de barras de tubos Fluent ID, verifique o LED de estado do leitor. Consulte a secção "LED de estado do Fluent ID" [> 93].
- Verifique a função do mostrador, a função do botão ou a mensagem de erro. Consulte as secções "Área de trabalho" [▶ 87] e "Botões de recuperação do método" [▶ 88].



- 4. Verifique a tabela da resolução de problemas. Consulte a secção "Tabelas da resolução de problemas" [▶ 166].
- 5. Se não for possível resolver o problema, consulte "Apoio ao cliente" [▶ 213].

6.6 Operação DeckCheck

Se o seu guião incluir a utilização de DeckCheck, o sistema de câmara DeckCheck irá captar imagens da mesa de trabalho depois do carregamento e comparar a disposição ao vivo com a disposição de referência.

O DeckCheck demora aproximadamente 20 segundos para um sistema de 3 braços/3 câmaras e aproximadamente 12 segundos para um sistema de 1 ou 2 braços/1 câmara para captar imagens da mesa de trabalho e apresentar a comparação das disposições ao vivo e de referência (assumindo que a configuração do PC é adequada – consulte o Manual de software de aplicação do FluentControl).

Tenha em atenção que na primeira utilização após ligar o instrumento, o comando DeckCheck irá precisar de mais tempo para apresentar o primeiro resultado – isto pode demorar alguns minutos.

Durante este tempo, o LED traseiro estará ligado.

Em sistemas Fluent de 3 braços, o braço central precisa de se mover entre as posições esquerda e direita (em sistemas de 1 ou 2 braços, os braços esquerdo e direito estarão posicionados no lado extremo esquerdo e extremo direito, respetivamente). Para esta movimentação do braço, a porta tem de estar fechada. Se a imagem for captada enquanto a porta está aberta, uma câmara estará geralmente bloqueada pelo braço central ou por qualquer braço que tenha sido movido manualmente.

Durante o processo de DeckCheck, o ecrã tátil apresenta formas a moverem-se no ecrã e a **Recolher imagens** seguido de **A verificar**. Após 12–20 segundos, dependendo do tamanho e da configuração do instrumento, as imagens da mesa de trabalho serão apresentadas em modo alternado. A imagem apresentada será a da primeira câmara com uma discrepância observada, começando pela esquerda.

O ecrã DeckCheck apresenta uma discrepância relativa à imagem de referência. Aqui, a imagem de referência mostra que deveria estar presente uma placa e que a câmara central detetou a discrepância.





Se lhe for apresento o botão **Ignorar e continuar**, o processamento continuará com a mesa de trabalho em vigor. Selecione o botão **Ignorar e continuar** se tiver a certeza de que não existem mais diferenças em relação à mesa de trabalho exigida, antes de fechar a porta. Selecione **Verificar** se pretender captar novas imagens da plataforma – tenha em atenção que se a porta não estiver fechada num sistema de 3 braços, a imagem será captada, mas o braço central irá obstruir a câmara. Caso contrário, será executada uma verificação automaticamente quando a porta estiver fechada.



Tenha em atenção que se o guião incluir a opção **mostrar sempre**, o ecrã acima será exibido e nenhuma discrepância será realçada. Contudo, as imagens ao vivo e de referência irão alternar e podem existir pequenas diferenças que não são captadas pelo sistema, mas que são visíveis a olho nu – por exemplo, algumas diferenças de cor, tubos/pontas únicos em falta ou pequenas deslocações laterais. Consulte os limites indicados abaixo.

Se forem detetadas discrepâncias, estas serão realçadas.

Para corrigir diferenças:

- 1. Abra a porta e substitua ou corrija a posição dos itens realçados.
- 2. O DeckCheck irá trabalhar continuamente para comparar a situação ao vivo corrigida com a disposição de referência.
- Utilize os botões DeckCheck para visualizar as diferenças captadas por cada uma das câmaras ou para pausar a visualização e conservar a imagem de referência conforme necessário. Quando não forem detetadas mais diferenças, o botão verde "Continuar" irá aparecer.
- 4. Selecione **Continuar** para prosseguir com o método.





Se algumas diferenças restantes forem, de facto, consideradas aceitáveis (por ex., o número total de pontas pode ser variável no início do método ou os níveis de líquido podem variar significativamente no início do processamento), pode selecionar a opção **Ignorar e continuar**, se disponibilizada no guião pelo operador principal.

Algumas diferenças de disposição podem não ser realçadas pelo DeckCheck – por ex., os seguintes tabuleiros de pontas coloridos do FCA:

Diferença entre os tipos de adaptadores de cabeças do MCA:

- Amarelo/laranja
- Branco/laranja
- Cinzento/todas as cores

Tipos de pontas diferentes do MCA 96

Tubos em falta em calhas de tubo parcialmente carregadas

Cartucho 300 SBS

Microplacas com rotação de 180 graus

Forma do poço da microplaca (por ex., poço redondo versus de fundo achatado ou de PCR)

Placas em hotéis periféricos com cartuchos de 10 ml/25 ml como aplique

Algumas tampas transparentes

Muitas destas diferenças são, contudo, claramente visíveis na mudança entre as disposições ao vivo e de referência.

6.7 Method Recovery

O FluentControl oferece a opção da recuperação de erros - por ex.:

A execução do método anterior foi abortada ou teve um erro fatal: a opção de recuperação do método oferece a possibilidade de continuar a partir do ponto em que ocorreu a falha na execução anterior.



Caso um método seja abortado ou tenha um erro fatal, deve ser realizada uma manutenção diária. Consulte os "Cuidados do sistema diários" [> 130].

6.7.1 Mudar para o modo de recuperação do método

- ✓ O operador principal ativou a opção de recuperação do método no FluentControl.
- ✓ A execução do método anterior foi abortada.



1. Selecione Method Recovery (A).



6.7.2 Recuperar a execução de um método

- ✓ A secção "Mudar para o modo de recuperação do método" [▶ 122] foi realizada.
- 1. Selecione a opção para continuar para o ecrã seguinte (A).

O ecrã apresenta a última linha de guia executada (C) e a linha de guia na qual a falha ocorreu – o ponto de recuperação (B).

Δ Method Recovery \otimes	
E Reagent Distribution	
Sample Transfer	— C
Get Tips	FCA 1 FCA, 200ul SBS
Aspirate	FCA 1 100 (µl) Water Video Single 1x24 13x75mm Tube Runner[001] Q1 - X1
Dispense	FCA 1 100 [µ1] Water Video Single 96 Well Flat[001] A1 - H1
Drop Tips	FCA-1 Wash Station Thru Deck DiTi Waste_1
🗉 🏣 Incubate & Wash	
Read & Export Data	
	A ×
Discard	Worktable Cancel

- Selecione os botões necessários descritos na secção "Botões de recuperação do método" [▶ 88].
- 3. Certifique-se de que a disposição da mesa de trabalho física do Fluent corresponde à disposição da mesa de trabalho (D) apresentada no ecrã tátil.
- 4. Selecione Run Recovery (E).



O sistema irá iniciar.



6.8 Desligar o instrumento

Caso não esteja nenhum método em execução, o instrumento passa para o modo de standby. O instrumento não necessita de ser desligado da rede.

Para desligar o instrumento, proceda do seguinte modo:

- 1. Pare qualquer método e selecione o modo de standby no ecrã tátil.
- 2. Certifique-se de que o instrumento está no modo de standby. Consulte a secção "Sinais de erro e estado do instrumento" [▶ 91].
- 3. Coloque os braços robóticos numa área de movimentação livre.





4. Desligue a energia utilizando o interruptor de energia (A) na traseira da fonte de alimentação externa.



7 System Care

Este capítulo dá instruções relativamente a todas as tarefas de cuidados do sistema a realizar para manter o Fluent em boas condições de funcionamento.



Opere o Fluent apenas quando este estiver em boas condições de funcionamento. Respeite rigorosamente as instruções de cuidados do sistema, conforme descritas neste manual. Para garantir o desempenho ideal e a fiabilidade, realize as tarefas de manutenção e de limpeza regularmente.

Na eventualidade de problemas ou questões, consulte a secção "Apoio ao cliente" [> 213].



O braço embate com objetos na mesa de trabalho

Os braços do Fluent podem ser movidos manualmente. Certifique-se de que os movimentos manuais dos braços são efetuados com suavidade, mantendo o braço fixo. Não deixe os braços embaterem com objetos sólidos, incluindo o batente mecânico do alcance do braço

7.1 Descontaminação

A descontaminação, de acordo com os regulamentos laboratoriais, é necessária nas circunstâncias indicadas na secção "Declaração de descontaminação" [> 38].

▲ ATENÇÃO

Contaminação!

Os resíduos de substâncias no Fluent podem causar ferimentos pessoais e afetar a integridade do processo.

 Descontamine o Fluent e os respetivos componentes e acessórios antes de qualquer interação.

O método de descontaminação deve ser definido pelo operador principal, com base no tipo de contaminante e no grau de contaminação. Neste capítulo, são disponibilizadas orientações sobre a seleção dos agentes de descontaminação e os modos de aplicação.



Para obter informações sobre o tratamento de Vapor de peróxido de hidrogénio, consulte o Manual de referência. Consulte os Documentos de referência.



Resultados de medição incorretos do Frida Reader!

Se não for montada nenhuma inserção, o Frida Reader pode fornecer resultados de medição incorretos.

 Utilize o bujão cego vermelho caso a inserção seja retirada (por exemplo, para limpeza).

7.2 Agentes de limpeza

7.2.1 Especificações dos agentes de limpeza

São necessários agentes de limpeza especiais para os cuidados do sistema. Todos os agentes de limpeza recomendados foram cuidadosamente selecionados e testados.

AVISO

Eficácia e compatibilidade química reduzida!

Não existe qualquer garantia relativamente à eficácia dos agentes de limpeza e à compatibilidade química caso sejam utilizados agentes de limpeza que não os recomendados pela Tecan.

- Utilize apenas os agentes de limpeza recomendados pela Tecan.
- Os agentes de limpeza são definidos para cada utilização específica nas tabelas dos cuidados do sistema. Não utilize agentes de limpeza salvo quando especificado para utilização numa tarefa específica.

A tabela seguinte especifica os agentes de limpeza especificados para utilização, conforme descrito nas tabelas de cuidados do sistema e nas atividades de cuidados do sistema:

Tab. 34: Agentes de limpeza para utilização conforme descrito nas tabelas de cuidados do sistema

Agente	Especificação
Água desionizada	Água destilada ou desionizada
Álcool	70% etanol, 100% isopropanol (2- propanol)
Detergente suave	Liqui-Nox
Agente ativo para superfícies	Contrad 70, Contrad 90/Contrad 2000, Decon 90
Desinfetante	Bacillol plus, SporGon
Desinfetante para superfícies (para contaminação por ácido nucleico)	DNAzap



Agente	Especificação
Ácido suave	ácido sulfúrico 0,3 M, ácido acético 10%, ácido fórmico 30-40%
Base	hidróxido de sódio 0,1 M
Lixívia	2% hipoclorito de sódio
Líquido do sistema	Conforme definido no método. Observe que as soluções aquosas com teor de sal devem ser purgadas durante a inatividade do sistema, por exemplo, durante a noite ou nos fins de semana. Consulte cuidados do sistema Final do dia.

7.2.2 Agentes de limpeza comerciais

Todas as instruções – indicadas pelo fabricante dos agentes de limpeza ou indicadas neste manual – para o manuseamento dos agentes de limpeza devem ser lidas atentamente e seguidas.

A tabela abaixo contém uma série de agentes de limpeza e desinfetantes disponíveis no mercado, especificados para utilização conforme descrito nas tabelas de cuidados do sistema e nas atividades de cuidados do sistema.

Agente de limpeza	Categoria do agente	Fabricante
DNAzap	Desinfetante para superfícies (para superfícies contaminadas com ácidos nucleicos)	Ambion www.ambion.com
Decon, Contrad	Agente ativo para superfícies	Decon Laboratories www.deconlabs.com
SporGon	Desinfetante	Decon Laboratories www.deconlabs.com
Bacillol Plus	Desinfetante	www.bode-chemie.com
Liqui-Nox	Detergente suave	Alconox www.alconox.com

Tab. 35: Agentes de limpeza comerciais

7.3 Modo de Cuidados do sistema

O operador principal define os métodos de cuidados do sistema necessários, de acordo com as tabelas de cuidados do sistema apresentadas em "Tabelas de cuidados do sistema" [▶ 130]. O modo de **System Care** faculta orientações para as tarefas de cuidados do sistema. Aceda a este modo através do ecrã tátil.



7.3.1 Mudar para o modo de Cuidados do sistema

- ✓ Os métodos de cuidados do sistema têm de estar disponíveis.
- 1. Selecione **System Care** (A).



2. Selecione a tarefa a executar.



- 3. Prima **Run** para iniciar o método de cuidados do sistema.
- 4. Realize as tarefas de cuidados do sistema.

7.3.2 Repor os erros

Se for apresentada uma mensagem, proceda do seguinte modo:

Mensagem

- Verifique a função do mostrador, a função do botão ou a mensagem de erro. Consulte as secções "Área de trabalho" [▶ 87] e "Botões de recuperação do método" [▶ 88].
- 2. Siga as instruções deste manual e do ecrã tátil para corrigir o erro.



 Continue a execução do método. Consulte a secção "Botões do mostrador, de opção e de ação" [▶ 88].

Se a lâmpada de estado acender ou mudar de cor, proceda do seguinte modo:

- 1. Verifique o estado do instrumento. Consulte a secção "Sinais de erro e estado do instrumento" [▶ 91].
 - Se o Fluent estiver equipado com um leitor de códigos de barras de tubos Fluent ID, verifique o LED de estado do leitor. Consulte a secção "LED de estado do Fluent ID" [▶ 93].
 - Verifique a função do mostrador, a função do botão ou a mensagem de erro. Consulte as secções "Área de trabalho" [▶ 87] e "Botões de recuperação do método" [▶ 88].
 - 4. Verifique a tabela da resolução de problemas. Consulte a secção "Tabelas da resolução de problemas" [▶ 166].
 - 5. Se não for possível resolver o problema, consulte "Apoio ao cliente" [▶ 213].

7.4 Tabelas de cuidados do sistema

Para garantir o desempenho ideal e a fiabilidade, realize as tarefas de manutenção e de limpeza conforme recomendado.



As tarefas das tabelas de cuidados do sistema apenas podem ser realizadas no modo de Cuidados do sistema. Consulte a secção "Modo de Cuidados do sistema" [> 128].

As tarefas de cuidados do sistema devem ser realizadas em intervalos regulares – nomeadamente, os cuidados do sistema diários, semanais e mensais.

7.4.1 Cuidados do sistema diários

7.4.1.1 Início do dia

Execute o método **Cuidados do sistema diários**, caso tenha sido disponibilizado pelo operador principal. Em alternativa, realize todas as tarefas individuais aplicáveis à configuração do seu braço do Fluent, apresentadas na tabela seguinte por ordem cronológica.

Tab. 36: Tabela de cuidados do sistema no início do dia

Instrumento/ Componente	Tarefa de cuidados do sistema	Agente de limpeza/ Produto descartável/ Dispositivo	Referência/ Atividades de cuidados do sistema
Pontas de perfuração	Inspecionar visualmente quanto à presença de acumulações nas pontas de perfuração. Limpar, se necessário. Verificar se as pontas não estão dobradas.	70% etanol ou 2% lixívia e pano sem fios	Consulte a secção "Limpar as pontas de perfuração" [▶ 143].

Lâmpada de estado



Instrumento/ Componente	Tarefa de cuidados do sistema	Agente de limpeza/ Produto descartável/ Dispositivo	Referência/ Atividades de cuidados do sistema
Cones de ponta descartável e pontas fixas	Inspecionar em relação a danos e acumulações	_	Esta tarefa está incluída no método Cuidados do sistema diários. AVISO! As pontas descartáveis não se destinam à reutilização.
Pontas fixas	Limpar. Verificar se as pontas não estão dobradas. Inspecionar visualmente com um espelho de dentista, para garantir que o revestimento está intacto.	70% etanol ou 100% isopropanol e pano sem fios	Consulte a secção "Limpar as pontas fixas" [▶ 143].
Recipiente de líquido do sistema (Liquid FCA com pontas fixas)	Garantir que está limpo e cheio, sem bolhas visíveis Certificar-se de que as tubagens para os conetores do recipiente estão unidas de forma adequada	_	Esta tarefa está incluída no método Cuidados do sistema diários.
Recipiente para resíduos do líquido (Liquid FCA com pontas fixas)	Garantir que está vazio Certificar-se de que as tubagens para os conetores do recipiente são novamente unidas de forma adequada	-	Esta tarefa está incluída no método Cuidados do sistema diários.
Saco de resíduos de pontas descartáveis	Garantir que está vazio	-	Consulte a secção "Substituir o saco de resíduos de pontas descartáveis" [▶ 155]. Esta tarefa está incluída no método Cuidados do sistema diários.



Instrumento/ Componente	Tarefa de cuidados do sistema	Agente de limpeza/ Produto descartável/ Dispositivo	Referência/ Atividades de cuidados do sistema
Sistema de líquido (Liquid FCA)	Garantir que está limpo	Líquido do sistema, álcool, água desionizada	Esta tarefa está incluída no método Cuidados do sistema diários . Em alternativa, pode ser executada separadamente como método Manutenção de purga de rotina do Liquid FCA . Consulte a secção "Limpar o caminho do líquido" [▶ 157].
Sistema de líquido (Liquid FCA)	Verificar visualmente a ausência de gotículas nas pontas ou cones DiTi após a purga	-	Esta tarefa está incluída no método Cuidados do sistema diários.
Dedos do manipulador	Verificar se os dedos estão direitos e nivelados Inspecionar em relação a danos e ao alinhamento incorreto	_	Em caso de desalinhamento, consulte a secção "Resolução de problemas do Robotic Gripper Arm (RGA)" [▶ 176]. Deformação ou danos. Consulte a secção "Apoio ao cliente" [▶ 213].
Dedos do manipulador FCA	Inspecionar em relação a danos	_	Substituir em caso de danos. Para obter informações de encomenda, consulte o manual de referência. Consulte a secção Documentos de referência.
Leitor Frida	Remova o bujão cego e encaixe a inserção no leitor Frida	-	Consulte a secção "Frida Reader" [▶ 162].

Resultados de medição incorretos do Frida Reader!

Se não for montada nenhuma inserção, o Frida Reader pode fornecer resultados de medição incorretos.

 Utilize o bujão cego vermelho caso a inserção seja retirada (por exemplo, para limpeza).

7.4.1.2 Final do dia

A tabela seguinte apresenta as tarefas de cuidados do sistema diárias do final do dia, por ordem cronológica:



Instrumento/ Componente	Tarefa de cuidados do sistema	Agente de limpeza/ Produto descartável/ Dispositivo	Referência/ Atividades de cuidados do sistema
Tabuleiros de recolha	Verificar se existem derramamentos e limpar ou substituir conforme necessário.	Água desionizada, álcool, detergente suave, desinfetante, base, lixívia, DNAzap	Consulte a secção "Limpar os tabuleiros de recolha da mesa de trabalho" [▶ 148].
Segmentos Caixa do Fluent ID	Limpar	Água desionizada, álcool, detergente suave, desinfetante, base, lixívia, DNAzap AVISO! Os agentes de limpeza necessários para a janela do leitor são diferentes dos necessários para o segmento em si. Consulte o "Cuidados do sistema semanais" [▶ 135].	Consulte a secção "Limpar as calhas e os segmentos" [▶ 148]. ATENÇÃO! Não olhe diretamente para o raio laser.
Película refletora (Fluent ID, Rotador de tubos)	Limpar e inspecionar em relação a danos	Álcool AVISO! Os agentes de limpeza necessários para a película refletora são diferentes dos necessários para o segmento em si.	Danos. Consulte a secção "Substituir a película refletora do Fluent ID" [▶ 150].
Calhas	Limpar	Água desionizada, álcool, detergente suave, desinfetante, agente ativo para superfícies, ácido suave, base, lixívia, DNAzap	Consulte a secção "Limpar as calhas e os segmentos" [▶ 148].
Pontas fixas	Limpar	Álcool, lixívia, pano sem fios	Consulte a secção "Limpar as pontas fixas" [▶ 143].
Cones de ponta descartável	Limpar	Álcool, pano sem fios	Consulte a secção "Limpar o cone de ponta descartável" [▶ 142].
Estação de lavagem e de resíduos (Liquid FCA)	Limpar	Água desionizada, álcool, detergente suave, desinfetante	Consulte a secção "Limpar os resíduos de pontas descartáveis e a unidade da estação de lavagem" [▶ 152].

Tab. 37: Tabela de cuidados do sistema no final do dia



Instrumento/ Componente	Tarefa de cuidados do sistema	Agente de limpeza/ Produto descartável/ Dispositivo	Referência/ Atividades de cuidados do sistema
Calha para resíduos de pontas descartáveis e tampas de resíduos	Limpar	Água desionizada, álcool, detergente suave, desinfetante	Consulte a secção "Limpar a calha para resíduos de pontas descartáveis" [▶ 153].
Sistema de líquido (Liquid FCA)	Purgar	Líquido do sistema AVISO! Se o sistema de líquido tiver um elevado teor de sal, purgue-o com água desionizada.	Execute o método Manutenção de purga de rotina do Liquid FCA.
Saco de resíduos de pontas descartáveis	Substituir	Especificações recomendadas do saco: L x C: 300 mm x 600 mm; espessura: 0,5 mm Material: polipropileno, polietileno ou co- polímero (autoclavável) AVISO! O saco de resíduos utilizado tem de cumprir as diretivas locais de segurança.	Consulte a secção "Limpar a calha para resíduos de pontas descartáveis" [▶ 153].
Recipiente de líquido do sistema (Liquid FCA)	Garantir que está limpo	Líquido do sistema	Consulte a secção "Conectar o recipiente do líquido do sistema e o recipiente para resíduos" [> 158].
Recipiente para resíduos (Liquid FCA com pontas fixas)	Esvaziar e limpar	Água desionizada, álcool, detergente suave, agente ativo para superfícies, desinfetante, base, lixívia	Dependendo das suas regras/regulamentos laboratoriais locais, limpe diária ou semanalmente. Consulte a secção "Conectar o recipiente do líquido do sistema e o recipiente para resíduos" [▶ 158].
Painel de segurança	Limpar	Água desionizada, álcool, detergente suave	Consulte a secção "Limpar os painéis de segurança" [▶ 152].
Rotador de tubos	Limpar as superfícies, o fixador e a estação de lavagem	Panos sem fios com 2% lixívia, 70% etanol ou 100% isopropanol	Consulte a secção "Limpar o Rotador de tubos" [▶ 143].



Instrumento/ Componente	Tarefa de cuidados do sistema	Agente de limpeza/ Produto descartável/ Dispositivo	Referência/ Atividades de cuidados do sistema
Pontas de perfuração	Inspecionar visualmente quanto à presença de acumulações nas pontas de perfuração. Limpar, se necessário.	70% etanol ou 2% lixívia e pano sem fios	Consulte a secção "Limpar pontas de perfuração" [▶ 143].
Estação de lavagem com ponta de perfuração e tubos de resíduos	Limpe o material de amostra residual da estação de lavagem e dos tubos.	Água desionizada, detergente suave, desinfetante. Não utilize lixívia sem enxaguar posteriormente os componentes da estação de lavagem com água.	Execute uma lavagem prolongada da estação de lavagem, incluindo todos os compartimentos (frontal, central e traseiro da estação de lavagem).
Leitor Frida	Remova a inserção e encaixe o bujão cego no leitor Frida	-	Consulte a secção "Frida Reader" [▶ 162].

7.4.2 Cuidados do sistema semanais

Os cuidados do sistema semanais devem ser realizados no último dia de trabalho de cada semana.

Execute o método **Cuidados do sistema semanais**, caso tenha sido disponibilizado pelo operador principal. Em alternativa, realize, além das tarefas diárias, todas as tarefas individuais aplicáveis à configuração do seu braço do Fluent, apresentadas na tabela seguinte por ordem cronológica.

Tab. 38: Tabela dos cuidados do sistema semanais

Instrumento/ Componente	Tarefa de cuidados do sistema	Agente de limpeza/ Produto descartável/ Dispositivo	Referência/ Atividades de cuidados do sistema
Sistema de líquido (Liquid FCA)	Limpar	Dependendo do líquido tratado pelo Fluent Decon, Contrad, base, ácido fraco, desinfetante	Consulte a secção "Limpar o caminho do líquido" [▶ 157].
		Seguido de purgas de água, álcool e líquido do sistema	



Instrumento/ Componente	Tarefa de cuidados do sistema	Agente de limpeza/ Produto descartável/ Dispositivo	Referência/ Atividades de cuidados do sistema
Liquid FCA	Verificar o aperto correto das seringas na interface da válvula e o aperto correto do êmbolo da seringa no parafuso de fixação do êmbolo.		Consulte a secção "Verificar o aperto das seringas" [▶ 159]
Cones de DiTi	Verificar o aperto do cone de DiTi	_	Consulte a secção "Apertar um cone DiTi" [▶ 161]
Liquid FCA	Realizar um teste de fugas (Liquid FCA)	_	Executar o Teste de fugas do Liquid FCA.
Air FCA	Realizar um teste de fugas (Air FCA)	_	Executar o Autoteste de cLLD e o teste de fugas do AirFCA.
Air FCA MultiSense	Executar o autoteste da cLLD do MultiSense Air FCA	_	Executar o Autoteste de cLLD e o teste de fugas do AirFCA.
Recipiente de líquido do sistema	Limpar	Água desionizada, álcool, detergente suave, agente ativo para superfícies, desinfetante, base, lixívia	Consulte a secção "Limpar o recipiente do líquido do sistema e o recipiente para resíduos" [▶ 159].
Estação de lavagem (Liquid FCA)	Limpar	Detergente ou solução antissética	-
Pontas dos dedos do manipulador RGA	Remover partículas e resíduos das pontas dos dedos do manipulador	Pano sem fios com álcool	Limpar com agente de limpeza.
Estação de encaixe e dedos do manipulador (interface de ligação)	Remover partículas e resíduos da interface de ligação do dedo do manipulador (PCBA, íman e cone)	Pano sem fios com álcool	Limpar com agente de limpeza.



Instrumento/ Componente	Tarefa de cuidados do sistema	Agente de limpeza/ Produto descartável/ Dispositivo	Referência/ Atividades de cuidados do sistema
Janela do leitor de código de barras autónomo	Limpar	Detergente suave	ATENÇÃO! Não olhe diretamente para o raio laser. Consultar o manual do fabricante do leitor de código de barras. Consultar a secção "Instrumento com radiação
Fluent ID e janela do leitor do rotador de tubos	Verificar se existe sujidade e danos Limpar, se necessário	Detergente suave Água desionizada para enxaguar	laser" [▶ 37]. ATENÇÃO! Não olhe diretamente para o raio laser. Limpar e enxaguar com um pano suave.
Fluent ID e refletor do rotador de tubos	Verificar se existe sujidade e danos Limpar, se necessário	Detergente suave Água desionizada para enxaguar	ATENÇÃO! Não olhe diretamente para o raio laser. Limpar e enxaguar com um pano suave.
Manipulador FCA	Limpar	Álcool	-
Rotador de tubos	Verificar a presença e o aperto dos pinos de bloqueio e de posicionamento. Apertar ou substituir os pinos, se necessário	-	Consultar a secção "Substituir pinos de bloqueio e pinos de posicionamento" [▶ 208].
MCA 96	Realizar um teste de fugas	-	Executar o método de fugas do MCA 96
MCA 96	Executar um teste de desempenho de pipetagem		 Executar o método de desempenho de pipetagem do MCA 96 se: A vida útil do vedante do cone tiver atingido os 90% Forem utilizadas pontas de 10 µl ou de 50 µl O Fluent Control emite um aviso para este efeito assim que forem atingidos 90% da vida útil do vedante do cone.



7.4.3 Cuidados do sistema mensais

A tabela seguinte apresenta as tarefas de cuidados do sistema mensais, por ordem cronológica:

Tab. 39: Tabela dos cuidados do sistema mensais

Instrumento/ Componente	Tarefa de cuidados do sistema	Agente de limpeza/ Produto descartável/ Dispositivo	Referência/ Atividades de cuidados do sistema
Software	Reiniciar o computador	_	Desligue o computador. Aguarde 10 segundos. Ligue o computador novamente.
Guia do braço	Limpar	Cotonete ou pano sem fios numa chave de fendas	Consulte a secção "Limpar a guia do braço" [▶ 160].
MCA 96	Limpar o MCH 96	Pano sem fios com álcool e ar comprimido	Consultar a secção "Limpar o MCH 96" [▶ 163]

7.4.4 Cuidados do sistema periódicos



Os intervalos para a realização destas tarefas devem ser determinados pelo operador principal.

A tabela seguinte apresenta as tarefas de cuidados do sistema, por ordem cronológica:

Tab. 40: Tabela dos cuidados do sistema periódicos

Instrumento/ Componente	Tarefa de cuidados do sistema	Agente de limpeza/ Produto descartável/ Dispositivo	Referência/ Atividades de cuidados do sistema
Conexão entre o cone e a manga	Remover partículas Limpar as superfícies	Álcool, pano sem fios	
Luz UVC	Verificar se existem impressões digitais. Limpar, se necessário.	Álcool, pano sem fios	
Dedos do manipulador do MCA 96	Inspeção visual dos dedos do manipulador, especialmente após colisões que os envolvam	Verificar os dedos do manipulador quanto a sujidade e danos. Em caso de sujidade, limpá-los com álcool e um pano sem fios.	



7.4.5 Cuidados do sistema anuais

Os cuidados do sistema anuais ajudam a manter a exatidão e precisão do instrumento, para minimizar os seus tempos de inatividade. Também ajudam a prolongar o período de vida do Fluent.

Contacte o responsável local pela assistência técnica da Tecan para agendar a marcação para os cuidados do sistema anuais. Consulte a secção "Apoio ao cliente" [> 213].

7.4.6 Cuidados do sistema bienais

As seguintes tarefas de manutenção preventiva têm de ser realizadas a cada 2 anos:

Tab. 41: Cuidados do sistema bienais

Componente	Tarefa	Referência
Manipulador FCA	Substituir dedos do manipulador FCA. Contador de reposições no FluentControl.	Para obter informações de encomenda, consulte o manual de referência. Consulte o "Documentos de referência" [▶ 9].

7.5 Atividades de Cuidados do sistema

Para realizar as atividades de cuidados do sistema descritas abaixo, proceda do seguinte modo:

- Mude para o modo de Cuidados do sistema. Consulte a secção "Modo de Cuidados do sistema" [▶ 128].
- Siga as instruções conforme descritas abaixo.

7.5.1 Deslocar o instrumento num armário dentro do laboratório

Danos no armário!

É possível que as prateleiras do armário tenham sido removidas, por exemplo, para a instalação da centrifugadora ou de um carrinho de resíduos. Deslocar o instrumento colocado num armário sem as respetivas prateleiras instaladas poderá danificar o armário e causar lesões.

- Antes de deslocar o instrumento, instale as prateleiras do armário.
- Desloque o armário apenas sobre um piso plano sem degraus ou sulcos. Na presença de degraus ou sulcos, utilize as barras de elevação do Fluent para elevar o sistema sobre a obstrução ou contacte o seu representante de assistência.

Para deslocar o instrumento num armário dentro de uma sala, proceda do seguinte modo:

1. Certifique-se de que o armário está estacionado em segurança e que está protegido contra o deslizamento.





2. Certifique-se de que as prateleiras do armário (A) se encontram instaladas.

- 3. Rode a porca do pé do armário (B) com uma chave de bocas.
- 4. Rode o parafuso vermelho (D) do pé do armário (C) até soltar o bloqueio e as rodas ficarem na posição móvel.



- 5. Mova o instrumento no armário até à nova localização.
- 6. Certifique-se de que o armário está estacionado em segurança e que está protegido contra o deslizamento.

7.5.1.1 Nivelar o instrumento

Para nivelar o instrumento, proceda do seguinte modo:





1. Com uma chave de bocas, baixe todos os pés ajustáveis até as rodas do

armário poderem ser rodadas manualmente.

2. Desaperte a porca de aperto (A) no pé correspondente.



3. Coloque o segmento de referência de acordo com as posições da grelha apresentadas abaixo.

Instrumento do tamanho 480: Posição da grelha do lado esquerdo 1 e posição da grelha do lado direito 21.

Instrumento do tamanho 780: Posição da grelha do lado esquerdo 1 e posição da grelha do lado direito 41.



Instrumento do tamanho 1080: Posição da grelha do lado esquerdo 1 e posição da grelha do lado direito 59.



- 4. Utilize o nível de bolha para garantir que o instrumento está alinhado horizontal e verticalmente.
- 5. Ajuste o nível do armário, conforme necessário (para a direita para levantar, para a esquerda para baixar).



- 6. Depois de nivelar o instrumento, volte a apertar as porcas de aperto nos pés do armário.
- 7. Certifique-se de que o armário está estacionado em segurança e que está protegido contra o deslizamento.

7.5.2 Limpar o cone de ponta descartável

Para limpar o cone de ponta descartável, proceda do seguinte modo:

- 1. Limpe os cones de pontas descartáveis com álcool, utilizando um pano sem fios.
- 2. Verifique os cones de pontas descartáveis e a ponta saliente durante os cuidados do sistema.

Para o Liquid FCA: Certifique-se de que a extensão da tubagem que sobressai do cone não está danificada.



3. Certifique-se de que as extensões da tubagem estão limpas e sem acumulações.

7.5.3 Limpar as pontas fixas

Riscos de lesões causadas pelas pontas fixas durante a limpeza

As pontas de pipetagem fixas podem provocar lesões.

 Quando estiver a utilizar a mesa de trabalho, evite o contacto com as pontas de pipetagem e com aerossóis, usando vestuário de proteção adequado.

Para limpar as pontas fixas, proceda do seguinte modo:

- 1. Limpe as pontas fixas com álcool, utilizando um pano sem fios.
- 2. Certifique-se de que as pontas fixas estão limpas e sem acumulações.

7.5.4 Limpar pontas de perfuração

Para limpar pontas de perfuração, realize o **método de Manutenção de limpeza de pontas de perfuração**. Este método tem de ser ajustado de acordo com a sua configuração da mesa de trabalho.

O guia inclui os seguintes passos:

- 1. Preparar a mesa de trabalho (isto é, material laboratorial e hardware).
- 2. Perfurar o Início Z de 8 tubos vazios com tampa num Rotador de tubos ou num carregador do fixador do tubo.
- 3. Limpar manualmente a parte acessível das pontas de perfuração com 70% etanol ou 2% lixívia com recurso a um pano sem fios. Evitar o contacto com o ápice afiado das pontas de perfuração.
- 4. Realizar comandos de lavagem após a limpeza manual.

7.5.5 Limpar o Rotador de tubos

Procedimento de limpeza geral

- 1. Para limpar qualquer parte do Rotador de tubos, utilize panos sem fios e impregne-os com um dos seguintes líquidos de limpeza: 2% lixívia, 70% etanol, 100% isopropanol
- Limpe as peças com panos embebidos para limpar e desinfetar. Utilize cotonetes para limpar áreas que não consegue alcançar com um pano sem fios.
- 3. Limpe os líquidos de limpeza com panos embebidos com água durante 5 minutos após a aplicação dos líquidos de limpeza.



Remover e limpar a placa do fixador do tubo

1. Para libertar a placa do fixador (A), segure na mesma com uma mão e puxe o pino de bloqueio do fixador (B) com a outra mão.

2. Remova a placa do fixador do rotador de tubos.



3. Limpe a placa do fixador de acordo com as instruções gerais acima ou, em alternativa, o fixador do tubo pode ser incubado num banho com 2% lixívia por um período máximo de 2 horas.

Limpar as superfícies do Rotador de tubos

1. Limpe as superfícies acessíveis do Rotador de tubos de acordo com as instruções gerais acima.




2. Para alterar a posição do tambor manualmente, segure no tambor com uma mão e prima o botão de libertação magnética.

- 3. Rode o tambor manualmente e solte o botão de libertação magnética.
- 4. Rode o tambor até este fixar no ponto magnético.
- 5. Limpe as superfícies anteriormente inacessíveis de acordo com as instruções gerais acima.

Montar a placa do fixador

- 1. Coloque a placa do fixador na parte superior do tambor do rotador de tubos.
- 2. Pressione verticalmente a placa do fixador do instrumento com uma mão e empurre o cursor preto para fixar a placa do fixador na sua posição.





7.5.6 Limpar a estação de lavagem do rotador de tubos

Procedimento de limpeza geral

- A estação de lavagem pode ser limpa na mesa de trabalho ou pode ser desmontada para limpeza.
- ✓ Utilize um escovilhão ao invés de panos para um melhor acesso.
- Para limpar qualquer uma das peças da estação de lavagem, utilize panos sem fios e impregne-os com um dos seguintes líquidos de limpeza: 2% lixívia, 70% etanol, 100% isopropanol
- 2. Limpe as peças com panos embebidos para limpar e desinfetar.
- 3. Limpe os líquidos de limpeza com panos embebidos com água durante 5 minutos após a aplicação dos líquidos de limpeza.

Desmontar a estação de lavagem do rotador de tubos

1. Empurre a alavanca de libertação da estação de lavagem na direção da caixa do leitor de código de barras e eleve a estação de lavagem com a outra mão.







2. Desconecte a tubagem de resíduos e coloque os conetores nos suportes da tubagem de resíduos.

Montar a estação de lavagem do rotador de tubos

1. Conecte os conetores da tubagem de resíduos.







 Monte a estação de lavagem nos eixos de orientação (A) e pressione-a verticalmente na direção da placa base.
 Verifique se a alavança de libertação ençaixa devidamente e fixa a estação

Verifique se a alavanca de libertação encaixa devidamente e fixa a estação de lavagem no respetivo lugar.



7.5.7 Limpar as calhas e os segmentos

Para limpar as calhas e os segmentos, proceda do seguinte modo:

AVISO

Falha na deteção de líquido (cLLD)!

Possível falha na deteção de líquido (cLLD) devido ao contacto comprometido entre a calha e o segmento da mesa de trabalho.

Certifique-se sempre de que as calhas e os segmentos estão limpos e secos.

1. Remova as calhas da mesa de trabalho do instrumento.

Os segmentos e os recetáculos estão a ser limpos no local.

2. Limpe a superfície das calhas, segmentos e recetáculos com um agente de limpeza.

Enxague as calhas, segmentos e recetáculos com água desionizada.

3. Volte a colocar as calhas na mesa de trabalho do instrumento.

7.5.8 Limpar os tabuleiros de recolha da mesa de trabalho

Para limpar os tabuleiros de recolha, proceda do seguinte modo:

- ✓ Os segmentos acima do tabuleiro de recolha são removidos. Consulte o "Remover segmentos" [▶ 99].
- ✓ Se não for possível remover segmentos da plataforma, como o Fluent ID, deslize os tabuleiros da plataforma para uma posição de plataforma aberta.
- 1. Remova os tabuleiros de recolha do instrumento.



- 2. Esvazie os tabuleiros, removendo o líquido de acordo com o protocolo laboratorial de manuseio do líquido para esse líquido.
- 3. Se os tabuleiros de recolha estiverem danificados ou em falta, devem ser substituídos.
- 4. Limpe a superfície dos tabuleiros de recolha com um agente de limpeza.
- Volte a colocar os tabuleiros de recolha no instrumento.
 Oriente os tabuleiros de recolha conforme ilustrado abaixo.
 Os tabuleiros de recolha adjacentes têm de ficar entrelaçados.



Fig. 41: Colocação incorreta do tabuleiro de recolha



Fig. 42: Tabuleiros de recolha entrelaçados





Fig. 43: Colocação correta do tabuleiro de recolha

7.5.9 Substituir a película refletora do Fluent ID

- ✓ Película refletora autocolante
- 1. Aqueça a película refletora. Utilize uma pistola de ar quente.
- 2. Remova a película refletora.



3. Remova os eventuais resíduos com álcool.





4. Aplique a nova película refletora autocolante na extremidade superior do refletor.

7.5.10 Aplicar a película refletora do Fluent ID na calha de resíduos DiTi

- ✓ Película refletora autocolante
- 1. Aplique fita na calha de resíduos DiTi, de acordo com a ilustração abaixo.



2. Aplique a nova película refletora na calha de resíduos DiTi, de acordo com a ilustração abaixo.





O raio laser deve estar no centro da película refletora.

3. Remova a fita da calha de resíduos DiTi.

7.5.11 Limpar os painéis de segurança

Para limpar os painéis de segurança, proceda do seguinte modo:

1. Limpe a superfície interior e exterior dos painéis de segurança com um agente de limpeza.

7.5.12 Limpar os resíduos de pontas descartáveis e a unidade da estação de lavagem

Para limpar os resíduos de pontas descartáveis e a unidade da estação de lavagem, proceda do seguinte modo:

- 1. Pressione o botão do fecho de libertação rápida (B).
- 2. Faça deslizar a estação de lavagem para trás.



Fig. 44: Fecho para a remoção da caixa do saco

3. Remova a estação de lavagem dos resíduos de pontas descartáveis e da unidade da estação de lavagem.





Fig. 45: Remoção da estação de lavagem

- 4. Limpe a superfície da estação de lavagem com um agente de limpeza e remova qualquer reagente derramado.
- 5. Pressione o botão do fecho de libertação rápida (B).
- 6. Coloque a estação de lavagem (A) na respetiva posição.
- 7. Empurre a estação de lavagem para a frente.



Fig. 46: A estação de lavagem é novamente colocada

7.5.13 Limpar a calha para resíduos de pontas descartáveis

Para limpar a calha para resíduos de pontas descartáveis, proceda do seguinte modo:

- ✓ O painel de segurança frontal está aberto.
- 1. Remova a tampa (A) da calha para resíduos de pontas descartáveis.





Fig. 47: Remoção da tampa da calha para resíduos de pontas descartáveis

2. Remova a calha para resíduos de pontas descartáveis (B) do suporte.



Fig. 48: Remoção da calha para resíduos de pontas descartáveis

3. Segure um tecido por baixo da abertura inferior da calha para resíduos de pontas descartáveis (C).

Impede o gotejar de substâncias contaminadas.



Fig. 49: Manuseio da calha para resíduos de pontas descartáveis

- 4. Coloque a calha para resíduos de pontas descartáveis e a tampa num recipiente cheio com agente de limpeza.
- 5. Deixe impregnar durante 30 minutos a 4 horas.
- 6. Retire a calha para resíduos de pontas descartáveis e a tampa do recipiente e coloque-os numa toalha limpa e seca.



- 7. Deixe secar.
- 8. Volte a colocar a calha para resíduos de pontas descartáveis (B) no suporte.



- Fig. 50: A inserção da calha para resíduos de pontas descartáveis é novamente colocada
- 9. Certifique-se de que o pino de posicionamento está corretamente colocado na ranhura (D).
- 10. Coloque a tampa (A) sobre a calha de resíduos.



Fig. 51: Pino de posicionamento e tampa

7.5.14 Substituir o saco de resíduos de pontas descartáveis

Para substituir o saco de desperdícios de pontas descartáveis, proceda do seguinte modo:

1. Levante o fecho (A) e faça deslizar a caixa do saco para a frente.







Fig. 52: Fecho para a remoção da caixa do saco

- 2. Remova a caixa do saco descartável (A).
- 3. Remova o saco de desperdícios de pontas descartáveis (B).
- 4. Elimine o saco de desperdícios de pontas descartáveis de acordo com as diretivas do seu laboratório.
- 5. Insira o saco de desperdícios descartável novo (B) na caixa do saco vazia (B).



Fig. 53: Caixa do saco e saco de desperdícios de pontas descartáveis





Fig. 54: Montagem correta da calha para resíduos na mesa de trabalho



Uma calha para resíduos mal posicionada pode provocar um embate com o braço e/ou a ejeção incorreta das pontas. Certifique-se de que as calhas para resíduos estão posicionadas corretamente, conforme ilustrado abaixo: corrija conforme o posicionamento da calha para resíduos

6. Faça deslizar a caixa do saco para a posição e vede-a com o fecho (A).



Fig. 55: Fecho de vedação

7.5.15 Limpar o caminho do líquido

- Para remover resíduos de proteínas do interior das pontas fixas, utilize ácido suave seguido de um agente de limpeza base.
- Para remover resíduos de ácido nucleico do interior das pontas fixas, utilize um agente de limpeza base.
- Agentes de limpeza, como o Decon/Contrad, podem afetar o processo. Portanto, caso sejam utilizados tais agentes, valide o processo cuidadosamente.
- Isopropanol é um desinfetante altamente eficaz. Evapora rapidamente, deixando as superfícies prontas para utilizar.
- Utilize apenas agentes de limpeza permitidos. Não utilize soluções de lixívia para purgar todo o sistema de líquido.

Para limpar o caminho do líquido, proceda do seguinte modo:

1. Separe a tubagem do líquido do sistema do recipiente de líquido.



- 2. Conecte o tubo de manutenção (30043739) à tubagem do sistema.
- 3. Insira a extremidade aberta da tubagem de manutenção numa garrafa com agente de limpeza.
- 4. Purgue com um agente de limpeza (20 ml com RapidWash e 10 ml com diluidor).
- 5. Deixe impregnar durante 20 minutos.
- 6. Coloque a tubagem numa garrafa com água desionizada.
- 7. Enxague duas vezes com água desionizada (20 ml com RapidWash e 10 ml com diluidor).

▲ ATENÇÃO

Líquidos inflamáveis!

Perigo de incêndio causado por líquidos inflamáveis ou líquido do sistema.

- Evite a formação e a acumulação de vapores inflamáveis.
- Não opere o sistema sem os tabuleiros de recolha.
- 8. Coloque a tubagem numa garrafa com álcool.
- 9. Purgue com álcool (20 ml com RapidWash e 10 ml com diluidor).
- 10. Remova a tubagem de manutenção da tubagem do sistema e conecte a tubagem do sistema ao recipiente do líquido do sistema.
- 11. Enxague duas vezes com água desionizada (20 ml com RapidWash e 5 vezes o volume do diluidor).
- 12. Verifique se existem bolhas na tubagem.
- 13. Purgue novamente se existirem bolhas.

7.5.16 Conectar o recipiente do líquido do sistema e o recipiente para resíduos

Para preparar o recipiente do líquido do sistema e o recipiente para resíduos, proceda do seguinte modo:



Só é garantido um funcionamento sem problemas se forem utilizados os recipientes originais com o sistema de controlo da Tecan. Antes da primeira utilização, o recipiente do líquido do sistema deve ser enxaguado manualmente e de forma meticulosa para remover quaisquer detritos sólidos do interior da garrafa. Consulte a secção "Limpar o recipiente do líquido do sistema e o recipiente para resíduos" [▶ 159].

- ✓ Recipiente da Tecan com uma capacidade de mais de 20 litros
- 1. Certifique-se de que o sistema de deteção do líquido (A, B) se encontra conectado corretamente.





2. Certifique-se de que os tubos (C, D) se encontram conectados corretamente.

7.5.17 Limpar o recipiente do líquido do sistema e o recipiente para resíduos

Para limpar o recipiente de líquido e o recipiente para resíduos, proceda do seguinte modo:

- 1. Esvazie o recipiente do líquido de lavagem manualmente.
- 2. Limpe o recipiente de líquido num reservatório com o agente de limpeza e enxague.
- 3. Desinfete o recipiente de líquido com álcool.
- 4. Conecte o líquido do sistema e o recipiente para resíduos. Consulte a secção "Conectar o recipiente do líquido do sistema e o recipiente para resíduos"
 [▶ 158].

7.5.18 Verificar o aperto das seringas

Para verificar o aperto correto das seringas, proceda do seguinte modo:





Fig. 56: Verificar o aperto

- A Parafuso de fixação do êmbolo B Parafuso da seringa
- Mova os êmbolos para o centro das seringas utilizando um script de manutenção que aspira ar.
 Nota: o script de manutenção tem de ser fornecido pelo administrador do FluentControl do laboratório.
- 2. Aperte o parafuso da seringa (B). Isto é, rode para a direita.
- 3. Rode o parafuso de fixação do êmbolo (A) para a direita para o apertar.

7.5.19 Limpar a guia do braço

Para limpar o a guia do braço, proceda do seguinte modo:

- 1. Limpe o rolamento da guia do braço (A) na respetiva guia com uma cotonete ou um pano sem fios numa chave de fendas.
- 2. Limpe os carris do braço (B) com um pano sem fios.





3. Quando aplicável, limpe a superfície superior do carril da guia da guia do braço MCA com um pano sem fios.

7.5.20 Apertar um cone DiTi

Para apertar o cone DiTi FCA, proceda do seguinte modo:

- 1. Segure o adaptador de pontas (D) e o tubo de ejeção de pontas (C).
- 2. Aperte o cone DiTi (A) com a chave para cones DiTi (B).



3. Execute o método de Manutenção de rotina do FCA .

7 - System Care Atividades de Cuidados do sistema



7.5.21 Frida Reader

Inserção



Fig. 57: Inserção do Frida Reader Para instalar, encaixe a inserção no Frida Reader e alinhe as marcas.

Bujão cego



Fig. 58: Bujão cego do Frida Reader

O bujão cego protege o Frida Reader quando a inserção é removida. Para instalar, encaixe o bujão cego no Frida Reader.



7.5.22 Limpar o MCH 96

Inspecione a placa de ejeção (A) para verificar a existência de danos visíveis. O derrame de determinados líquidos de processo (como DMSO ou acetonitrilo) ou agentes de limpeza (como lixívia) pode provocar danos na placa. Caso isto aconteça, contacte o serviço de assistência local para proceder à substituição.



- 1. Utilize a ferramenta de deslocação para deslocar o eixo de ejeção para a parte inferior. 2. . 3.
- 2. Utilize ar comprimido sem óleo para limpar a superfície superior da placa de ejeção (C), a placa de cone e os cones (B)
- Utilize um pano sem fios e álcool para limpar a superfície inferior da placa de ejeção (D).





A imagem em baixo à esquerda apresenta a placa de ejeção antes da limpeza e a imagem à direita apresenta a placa de ejeção após a limpeza.





8 Resolução de problemas

Consulte este capítulo para obter ajuda para retomar a operação depois de ter ocorrido um erro no Fluent. Para obter mais informações ou, em caso de problemas não abordados neste manual ou de detalhes insuficientes, consulte a secção "Apoio ao cliente" [> 213].

8.1 Instruções de segurança deste capítulo

Contaminação cruzada devido a pontas danificadas após embate!

As pontas dobradas ou o revestimento da ponta danificado causam imprecisões na pipetagem e erros na deteção do líquido.

Verifique as pontas fixas após um embate. Consulte a secção "Verificar as pontas fixas" [▶ 187].

8.2 Tabelas da resolução de problemas



As tabelas da resolução de problemas apresentam os possíveis problemas, causas e medidas corretivas. Para obter mais informações ou, em caso de problemas não abordados neste manual ou de detalhes insuficientes, consulte a secção "Apoio ao cliente" [> 213].

8.2.1 Resolução de problemas do instrumento

Problema/erro	Causa possível	Medida corretiva
Fuga de líquido do sistema	A tubagem e/ou as conexões da tubagem não estão apertadas. A seringa tem uma fuga.	Consulte a secção "Apoio ao cliente" [▶ 213].
Erro de comunicação	A energia não está ligada. A alimentação ou a comunicação foi interrompida. Sem comunicação.	Desligue o instrumento. Aguarde até a lâmpada de estado do instrumento e a lâmpada da fonte de alimentação desligarem. Desligue o PC. Verifique o cabo e as fichas. Ligue o instrumento e o PC.
	Acionamento X, Y ou Z bloqueado.	Verifique se existem obstáculos. AVISO! Certifique-se de que os braços podem mover-se livremente.

Tab. 42: Tabela da resolução de problemas do instrumento



Problema/erro	Causa possível	Medida corretiva
Erro de inicialização	Não é possível inicializar os braços.	Verifique se existem obstáculos. AVISO! Certifique-se de que os braços podem mover-se livremente.
	Problema de hardware.	Consulte a secção "Apoio ao cliente" [▶ 213].
O sensor da porta do painel de segurança frontal e o bloqueio da porta estão danificados	Falha mecânica dos bloqueios da porta.	Desligue o instrumento. Consulte a secção "Apoio ao cliente" [▶ 213].
Painel de segurança em falta ou danificado	Não é possível garantir a segurança.	Desligue o instrumento. Consulte a secção "Apoio ao cliente" [▶ 213].
Falha na deteção de líquido (cLLD)	Superfície de contacto suja. Contacto inadequado entre o material laboratorial e o segmento.	 Prepare a mesa de trabalho. Consulte a secção . Limpe a superfície de contacto. Consulte a secção "Limpar as calhas e os segmentos" [▶ 148].
	O líquido do sistema tem uma condutividade > 10 μS/cm para a compatibilidade da cLLD.	Contacte o operador principal.

8.2.2 Resolução de problemas do Flexible Channel Arm (FCA)

Problema/erro	Causa possível	Medida corretiva
Cone DiTi solto CUIDADO! Volu mes de pipetagem imprecisos!	Cone de DiTi apertado de forma insuficiente.	Apertar o cone de DiTi.
Ponta descartável não recolhida	Cone de DiTi apertado de forma insuficiente.	Apertar o cone de DiTi.

Tab. 43: Tabela da resolução de problemas do Flexible Channel Arm



Problema/erro	Causa possível	Medida corretiva
Ponta descartável não eliminada	Cone de DiTi apertado de forma insuficiente.	Apertar o cone de DiTi.
	DiTi reutilizadas	Garantir que as pontas são novas. Não é recomendada a reutilização de DiTi.
		Pontas descartáveis não ejetadas na calha para resíduos/a calha para resíduos não está posicionada corretamente
As pontas descartáveis não foram ejetadas para a calha para resíduos	A calha para resíduos não está posicionada corretamente	Certifique-se de que as calhas para resíduos estão posicionadas corretamente. Consulte a secção "Limpar a calha para resíduos de pontas descartáveis" [▶ 153]
Pontas não alinhadas com o material	Transportador na posição errada. O segmento não se	Garantir a posição correta do transportador. Consultar a secção "Carregar as calhas standard" [▶ 100].
laboratorial num transportador	encontra bloqueado no seu lugar. Material laboratorial posicionado incorretamente.	Bloquear o segmento no seu lugar. Consultar a secção "Verificar o segmento" [▶ 206].
Pontas não alinhadas com o material laboratorial em vários transportadores	Alinhamento incorreto do braço devido a uma colisão.	Consultar a secção "Apoio ao cliente" [▶ 213].
A ponta colide com a parte inferior do material laboratorial	Material laboratorial errado. Material laboratorial posicionado incorretamente.	Garantir que o material laboratorial na plataforma corresponde à disposição da plataforma do método.
A DiTi pinga	O cone de DiTi sujo causa fugas.	Limpar o cone de DiTi.
	DiTi reutilizadas	Garantir que as pontas são novas. Não é recomendada a reutilização de DiTi.
Mensagem de erro: Pressure out of range (Pressão fora do intervalo) (Air FCA)	Filtro em linha molhado após a aspiração com o tamanho de DiTi errado.	Garantir de que o tamanho da DiTi na plataforma corresponde ao definido para o método. Certifique-se de que os cones de DiTi estão apertados corretamente Verificar o filtro em linha. Consultar a secção "Verificar o filtro interno (Air FCA)" [▶ 180].



Problema/erro	Causa possível	Medida corretiva
Mensagens de erro:	Campo magnético a interferir com o sensor	-
DiTi not fetched (DiTi não recolhida)	de presença de DiTi.	
DiTi not dropped (DiTi não baixada)		
Deteção do líquido: A ponta não deteta líquido	Cone de DiTi solto	Apertar o cone de DiTi (consultar a secção "Apertar um cone DiTi" [▶ 161]).
O-rings gastos	Adaptadores de pontas MultiSense	Substituir os O-rings e o X-ring nos adaptadores de pontas MultiSense.

8.2.3 Mistura e perfuração

Tab. 44: Resolução de problemas

Sintoma	Causa possível	Medidas corretivas
A ponta de perfuração não pode ser recolhida através de comandos de software	Ponta de perfuração presa	Consulte a secção "Recolher pontas de perfuração presas" [▶ 198].
Ponta de perfuração danificada	Ponta de perfuração dobrada Ponta danificada	Substitua a ponta de perfuração. Consulte as secções "Remover pontas de perfuração" [▶ 191] e "Instalar pontas de perfuração" [▶ 194].



Sintoma	Causa possível	Medidas corretivas
Erros de perfuração	Ponta de perfuração demasiado seca	Lubrificação com água (estação de lavagem)
	Ponta danificada	Substitua a ponta de
	Ponta de perfuração dobrada	perfuração. Consulte as secções "Remover pontas de perfuração" [▶ 191] e "Instalar pontas de perfuração" [▶ 194].
	Parâmetro de perfuração errado	Contacte o operador principal.
	Tipo de movimento utilizado errado	Contacte o operador principal.
	Tubos utilizados errados	Utilize tubos compatíveis. Consulte a secção "Calhas do rotador de tubos" [▶ 78].
	O braço atingiu a sua vida útil	Consulte a secção "Apoio ao cliente" [▶ 213].
Problema com a manipulação de líquidos	Pontas de perfuração obstruídas	Lave as pontas de perfuração. Verifique o procedimento de lavagem na sua generalidade.
	Ponta danificada	Substitua a ponta de perfuração. Consulte as secções "Remover pontas de perfuração" [▶ 191] e "Instalar pontas de perfuração" [▶ 194].
	Seringas montadas de forma não adequada	Verifique o aperto das seringas. Consulte a secção "Verificar o aperto das seringas" [▶ 159].
	Bolhas no sistema de líquido	Purgar. Consulte a secção "Apoio ao cliente" [▶ 213].





Sintoma	Causa possível	Medidas corretivas
Problemas de hemólise	Diluição da amostra	Volume excessivo ou volume de repartição superior
		Solução salina de 0,9% como volume de repartição
		Velocidades de pipetagem menores
	Ponta danificada	Substitua a ponta de perfuração. Consulte as secções "Remover pontas de perfuração" [▶ 191] e "Instalar pontas de perfuração" [▶ 194].
	Parâmetros de mistura	Certifique-se de que os parâmetros de rotação/ oscilação utilizados do rotador de tubos não provocam hemólise



Sintoma	Causa possível	Medidas corretivas
Amostra na seringa	Qualquer	Limpe o sistema. Consulte a secção "Limpar o caminho do líquido" [▶ 157].
	Bolsa de ar errada	Valide o procedimento de lavagem.
	Seringas montadas de forma não adequada	Verifique o aperto das seringas. Consulte a secção "Verificar o aperto das seringas" [▶ 159].
		Execute o método de fugas do FCA.
		Bolsa de ar inicial maior.
		Velocidades de aspiração menores.
	Preparação incorreta das amostras para manipulação de líquidos. Os tubos de origem da amostra contêm partículas sólidas, como coágulos, detritos celulares, etc.	Garanta uma preparação de amostra adequada para permitir a pipetagem do líquido de amostra.
		Certifique-se de que os tubos de origem da amostra não contêm partículas sólidas, como coágulos, detritos celulares, etc.
	Preparação incorreta das amostras para manipulação de líquidos. Os tubos não estão devidamente cheios e ainda possuem um vácuo parcial que reduz a bolsa de ar inicial durante a perfuração.	Certifique-se de que os tubos de origem da amostra estão devidamente cheios com o volume alvo do tubo.
		Certifique-se de que os tubos de origem da amostra não contêm vácuo.
		Aumente a bolsa de ar inicial para compensar um possível vácuo residual.
A ponta de perfuração dobra-se durante o procedimento de lavagem	A ponta de perfuração não está centrada nos orifícios de limpeza da estação de lavagem.	Crie uma cópia da estação de lavagem e programe as posições de pipetagem.



Sintoma	Causa possível	Medidas corretivas
Deteção incorreta do nível de líquido: apenas em canais específicos	Ponta de perfuração dobrada: A ponta de perfuração está dobrada e, por isso, toca na parede do tubo durante a perfuração	Substitua a ponta de perfuração. Consulte as secções "Remover pontas de perfuração" [▶ 191] e "Instalar pontas de perfuração" [▶ 194].
	A posição da ponta está incorreta e, por isso, a ponta de perfuração toca na parede do tubo durante a perfuração.	Utilize material laboratorial fabricado pela Tecan. Consulte a secção "Calhas do rotador de tubos" [▶ 78].
		Programe/ajuste a posição de pipetagem do material laboratorial
	A orientação da ponta de perfuração está incorreta.	Monte a ponta de perfuração com a abertura virada para a frente do instrumento. Consulte a secção "Instalar pontas de perfuração" [▶ 194].
Deteção incorreta do nível de líquido: desvio constante do nível de líquido esperado e do nível de líquido detetado	As tolerâncias de fabrico do braço, do rotador de tubos e/ ou do instrumento em conjunto com a força de perfuração podem dar origem a um desvio Z significativo durante a deteção do nível de líquido.	Programe/ajuste o atributo personalizado "CompensaçãoAlturaDeteção Perfuração" na definição do material laboratorial do tubo



Sintoma	Causa possível	Medidas corretivas
Transbordo da estação de lavagem	Amostra obstruída na estação de lavagem, resíduos centrais, produtos de limpeza frontais ou posteriores	Remova a estação de lavagem e limpe conforme descrito. Consulte a secção "Limpar a estação de lavagem do rotador de tubos" [▶ 146].
		Implemente um procedimento de limpeza mais rigoroso como parte da manutenção no final do dia. Consulte as secções "Final do dia" [▶ 132] e "Mistura e perfuração" [▶ 169] (passo 6).
	Conetores de resíduos obstruídos	Limpe os conetores com a escova para garrafas ou substitua os conetores. Consulte a secção "Limpar a estação de lavagem do rotador de tubos" [▶ 146].
		Recomenda-se a substituição regular dos conetores da estação de lavagem. Consulte a secção "Instruções de segurança deste capítulo" [> 94] (aplicações de sangue total)
		Implemente um procedimento de limpeza mais rigoroso como parte da manutenção no final do dia. Consulte as secções "Final do dia" [▶ 132] e "Mistura e perfuração" [▶ 169] (passo 6).

8.2.4 Resolução de problemas do Multiple Channel Arm (MCA)

Tab. 45: Tabela da resolução de problemas do Multiple Channel Arm

Problema/erro	Causa possível	Medida corretiva
As pontas não estão alinhadas com os transportadores	Falha mecânica	Consulte a secção "Apoio ao cliente" [▶ 213].
	Falha do braço	
A microplaca e a cabeça de pipetagem não estão totalmente paralelas	Embate	Contacte o operador principal para verificar o paralelismo da cabeça de pipetagem e da plataforma.



Problema/erro	Causa possível	Medida corretiva
Durante a pipetagem, a cabeça de pipetagem para, gerando um erro	A aceleração de aspiração e de distribuição é demasiado rápida em comparação com a velocidade. A desaceleração de aspiração e de distribuição é demasiado rápida em comparação com a velocidade.	 A aceleração tem de ser proporcional à velocidade de aspiração e de distribuição. A desaceleração tem de ser proporcional à velocidade de aspiração e de distribuição. O problema não pode ser resolvido. Consulte a secção "Apoio ao cliente" [▶ 213].
Vários ou todos os canais de pipetagem com fugas	Pontas descartáveis ou vedações do cone de ponta erradas	Utilize sempre pontas descartáveis ou vedações do cone de ponta fornecidas pela Tecan. Contacte o operador principal para verificar a existência de fugas no sistema.
	As vedações do cone de ponta são antigas ou defeituosas.	Consulte a secção "Apoio ao cliente" [▶ 213]. Contacte o operador principal para verificar a existência de fugas no sistema.
	A cabeça de pipetagem está defeituosa.	Consulte a secção "Apoio ao cliente" [▶ 213].
Um único canal com fuga	A vedação do cone de ponta ou outras vedações da cabeça de pipetagem estão defeituosas.	Contacte o operador principal para verificar a existência de fugas no sistema. Consulte a secção "Apoio ao cliente" [▶ 213].
Uma única ponta descartável não recolhida corretamente	A ponta descartável individual está defeituosa. A vedação do cone de ponta nesta posição de ponta descartável está defeituosa.	Substitua as pontas descartáveis. O problema não pode ser resolvido. Consulte a secção "Apoio ao cliente" [▶ 213].
Uma ponta descartável não largada	A ponta descartável individual está defeituosa. A vedação do cone de ponta nesta posição de ponta descartável está defeituosa.	O problema não pode ser resolvido. Consulte a secção "Apoio ao cliente" [▶ 213].
	Humidade incorreta	Certifique-se de que a humidade se encontra dentro dos limites de humidade de funcionamento. Consulte a secção "Condições ambientais" [▶ 45].
Várias ou todas as pontas descartáveis ficaram por largar	Foram utilizadas pontas descartáveis erradas.	Utilize sempre as pontas descartáveis fornecidas pela Tecan. O problema não pode ser resolvido. Consulte a secção "Apoio ao cliente" [▶ 213].



Problema/erro	Causa possível	Medida corretiva
A caixa de pontas descartáveis é levantada com as pontas descartáveis ao recolher as pontas descartáveis	O transportador não está corretamente ajustado. O desvio X e/ou Y estão especificados incorretamente.	Ajuste todos os transportadores (mecânicos) com precisão. Substitua o transportador de pontas descartáveis. O problema não pode ser resolvido. Consulte a secção "Apoio ao cliente" [▶ 213].
	A caixa de pontas descartáveis não cumpre as especificações.	Utilize sempre caixas de pontas descartáveis que cumpram as normas da Society of Biomolecular Screening. O problema não pode ser resolvido. Consulte a secção "Apoio ao cliente" [▶ 213].
	O transportador de pontas descartáveis está defeituoso (falha nos retentores da caixa de pontas descartáveis).	Utilize sempre caixas de pontas descartáveis que cumpram as normas da Society of Biomolecular Screening. O problema não pode ser resolvido. Consulte a secção "Apoio ao cliente" [▶ 213].
Resultados de pipetagem imprecisos	As pontas descartáveis não são recolhidas adequadamente. Os parâmetros de manuseio do líquido estão incorretos. Os suportes não estão ajustados corretamente. A cabeça de pipetagem está defeituosa.	Contacte o operador principal para verificar o guião da aplicação e os transportadores. Contacte o operador principal para verificar os parâmetros ambientais e a altura de distribuição.

8.2.5 Resolução de problemas do Robotic Gripper Arm (RGA)

Tab. 46: Tabela da resolução de problemas do Robotic Gripper Arm

Problema/erro	Causa possível	Medida corretiva
Microplaca não recolhida	Sem microplacas no carregador. Os dedos do manipulador não conseguem recolher a microplaca.	Coloque microplacas no carregador. Configure a posição do manipulador. Limpe os dedos do manipulador do RGA.
Ruídos estranhos durante o movimento do braço	Os componentes estão danificados ou gastos.	Consulte a secção "Apoio ao cliente" [▶ 213].



Problema/erro	Causa possível	Medida corretiva
Dedos excêntricos do manipulador mal alinhados	Embate dos dedos sobressalentes. Parafusos dos dedos insuficientemente apertados.	Alinhe os dedos excêntricos do manipulador. Consulte a secção "Verificar o alinhamento do dedo do manipulador" [▶ 200].
		Utilize uma chave de fendas dinamométrica para apertar os parafusos para 3 Nm, conforme descrito na secção "Alinhamento básico dos dedos do manipulador para dedos do manipulador do FES" [> 201].

8.2.5.1 Resolução de problemas do Robotic Gripper Arm com eixo Z longo (RGA-Z)

Problema/erro	Causa possível	Medida corretiva
Microplaca não recolhida	Sem microplacas no carregador. Os dedos do manipulador não conseguem recolher a microplaca.	Coloque microplacas no carregador. Configure a posição do manipulador. Limpe os dedos do manipulador do RGA.
	Os dedos do manipulador estão escorregadios.	Limpe os dedos do manipulador do RGA.
Ruídos estranhos durante o movimento do braço	Os componentes estão danificados ou gastos.	Consulte a secção "Apoio ao cliente" [▶ 213].

Tab. 47: Tabela da resolução de problemas do Robotic Gripper Arm com eixo Z longo

8.2.6 Resolução de problemas do sistema de lavagem

Tab. 48: Tabela da resolução de problemas do sistema de lavagem

Problema/erro	Causa possível	Medida corretiva
Transbordo da estação de lavagem	A estação de lavagem está bloqueada por pontas descartáveis ou algas.	Limpe a estação de lavagem. Consulte a secção "Limpar o recipiente do líquido do sistema e o recipiente para resíduos" [▶ 159].
	A tubagem de desperdícios está dobrada.	Verifique se existem dobras na tubagem. Consulte a secção "Verificar a tubagem do recipiente do líquido do sistema e do recipiente para resíduos" [▶ 106].
	Os tubos de resíduos estão obstruídos ou danificados.	Verifique a tubagem de desperdícios. Substitua a bomba de desperdícios, se necessário. Consulte a secção "Apoio ao cliente" [▶ 213].





Para a resolução de problemas em sistemas de mistura e perfuração e de estações de lavagem, consulte "Mistura e perfuração" [▶ 169].

8.2.7 Resolução de problemas do Fluent ID

Tab. 49: Tabela da resolução de problemas do Fluent ID

Problema/erro	Causa possível	Medida corretiva
Código de barras não lido	Etiqueta do código de barras não virada para o leitor.	Descarregue a calha de tubos e vire os tubos de forma que as etiquetas dos códigos de barras fiquem viradas para a esquerda. Volte a carregar a calha de tubos no Fluent.
	Calha carregada demasiado rápido.	Descarregue a calha de tubos e volte a carregá-la lentamente.
	Má qualidade da etiqueta.	Introduza o código de barras manualmente ou comunique o problema ao operador principal.
	A janela do leitor está suja.	Limpe a janela do leitor. Consulte a secção .
	O refletor está sujo.	Limpe o refletor. Consulte a secção .
	O tipo de código de barras ou o comprimento do código de barras não está predefinido para o método.	Comunique o problema ao operador principal.
Presença do tubo não detetada	Posição da etiqueta do código de barras muito baixa no tubo.	Comunique o problema ao operador principal.

8.2.8 Resolução de problemas de software

Tab. 50: Tabela da resolução de problemas de software

Problema/erro	Causa possível	Medida corretiva
Ecrã do início de sessão do utilizador não apresentado quando previsto.	A gestão do utilizador não foi ativada no FluentControl.	Contacte o operador principal para ativar a gestão do utilizador.
O utilizador não consegue iniciar sessão.	A palavra-passe está incorreta ou a conta está bloqueada.	Contacte o operador principal para repor a palavra-passe ou a conta.
As ações de assistência não foram todas concluídas. Surgem advertências em cada arranque do FluentControl.	As ações de assistência previstas não estão todas marcadas como concluídas na configuração do instrumento.	Consulte a secção "Apoio ao cliente" [▶ 213].



Problema/erro	Causa possível	Medida corretiva
O ecrã tátil não reage ao toque.	Controlador de software não instalado.	Contacte o administrador do computador para instalar os controladores do CD de instalação e configurar o ecrã tátil.
	Interface do ecrã tátil configurada incorretamente.	Abra as definições do controlador do ecrã tátil e certifique-se de que o ecrã tátil está mapeado corretamente.
A interface de toque não é apresentada no ecrã tátil.	O ecrã tátil não estava ligado no momento do arranque do software.	Ligue o instrumento e reinicie o software ou verifique as definições da ferramenta de toque no sistema de configuração do FluentControl.
Erro no arranque do FluentControl.	O FluentControl (SystemSW.exe) já está em execução em segundo plano (Gestor de tarefas).	Abra o Gestor de tarefas, processe SystemSW.exe e reinicie o FluentControl. Ou reinicie o computador.
O FluentControl não comunica com os dispositivos de hardware conectados.	O FluentControl não está configurado corretamente para comunicar com os dispositivos de hardware.	Contacte a pessoa responsável pela configuração do sistema para ativar o estado I/O dos dispositivos de hardware.





8.3 Atividades de resolução de problemas

8.3.1 Posição do material laboratorial

Para se certificar de que o material laboratorial está posicionado corretamente no recetáculo para um acesso preciso do braço, proceda do seguinte modo:

1. Coloque o material laboratorial (A) no recetáculo (B).



 Pressione suavemente o material laboratorial contra o posicionador estático (C).



- 3. Faça deslizar o posicionador deslizante diagonal (E) no sentido do material laboratorial ou na direção oposta, para o ajustar com precisão.
- 4. Faça deslizar os posicionadores deslizantes vertical e horizontal (D, F) no sentido do material laboratorial ou na direção oposta, para o fixar.
- 5. Levante o material laboratorial do recetáculo.

Certifique-se de que não há fricção ao colocar ou remover a placa.

8.3.2 Verificar o filtro interno (Air FCA)

Está instalado em cada canal um sistema de controlo para proteger os canais de pipetagem do Air FCA contra a aspiração excessiva de líquido.


- O método Manutenção de rotina do Air FCA inclui uma verificação do filtro interno, em que são detetados os filtros molhados, danificados ou posicionados incorretamente e ainda os filtros em falta.
- Execute o método Manutenção de rotina do Air FCA para verificar o filtro interno no interior do cone DiTi de um canal de pipetagem do Air FCA.

Em caso de erro, o filtro interno deve ser substituído. Consulte a secção "Substituir o filtro interno (Air FCA)" [▶ 181].

8.3.3 Substituir o filtro interno (Air FCA)

Para substituir o filtro interno, proceda do seguinte modo:

- ✓ Cone de ponta descartável descontaminado.
- ✓ Cone DiTi removido do canal. Consulte a secção "Remover o cone DiTi (Air FCA)" [▶ 182].



- Perfure o filtro em linha (B) lateralmente com a ferramenta de remoção do filtro (C).
- 2. Retire o filtro em linha com a ferramenta de remoção do filtro. Observe que o filtro pode estar contaminado com os líquidos do processo.





3. Elimine o filtro interno.



- Limpe o cone DiTi (A) com álcool.
 O cone DiTi deve estar limpo antes de ser novamente colocado.
- 5. Coloque o novo filtro em linha sobre uma superfície limpa e plana.
- Pressione o filtro em linha para dentro do cone DiTi.
 O filtro em linha não deve sobressair do cone DiTi.
- 7. Verifique o filtro em linha de acordo com o método definido pelo seu operador principal.

8.3.4 Remover o cone DiTi (Air FCA)

Para remover o cone DiTi (Air FCA), proceda do seguinte modo:

- ✓ Chave para cones DiTi disponível.
- 1. Desligue o instrumento.
- 2. Abra o painel de segurança frontal.
- 3. Levante manualmente todas as rodas Z até à posição mais elevada.
- 4. Desloque todas as rodas Z no sentido frontal do instrumento.
- 5. Afaste o mais possível as rodas Z.
- 6. Segure o adaptador de pontas (D) e o tubo de ejeção de pontas (C).





7. Desenrosque o cone DiTi (A) com a chave para cones DiTi (B).

8. Puxe o cone DiTi para baixo com cuidado.

Em alguns casos, o tubo de ejeção de pontas (C) ou o cilindro adaptador (B) podem estar ainda conectados ao cone DiTi (A). Consulte a secção "Montar o tubo de ejeção DiTi (Air FCA)" [▶ 183].



8.3.5 Montar o tubo de ejeção DiTi (Air FCA)

Para montar o tubo de ejeção DiTi (Air FCA), proceda do seguinte modo:

- ✓ O tubo de ejeção DiTi foi removido, de acordo com as instruções.
- ✓ Chave para cones DiTi disponível.





1. Introduza a manga de vedação (B) no cilindro adaptador (A).

2. Enrosque o cone DiTi (C) no cilindro montado. Certifique-se de que o O-ring preto não está visível, conforme ilustrado na imagem abaixo.



3. Insira o tubo ejetor de pontas (D) com o lado sem ranhura no cilindro montado, conforme ilustrado abaixo.



8.3.6 Instalar o cone DiTi (Air FCA)

Para instalar o cone DiTi Air FCA, proceda do seguinte modo:

- ✓ O cone DiTi está totalmente montado: Consulte a secção "Montar o tubo de ejeção DiTi (Air FCA)" [▶ 183].
- ✓ Chave para cones DiTi disponível.
- 1. Introduza o cilindro adaptador no tubo de ejeção de pontas (C).
- 2. Segure o adaptador de pontas (D) e o tubo de ejeção de pontas (C).





3. Enrosque o cone DiTi (A) com a chave para cones DiTi (B).

4. Execute o método Manutenção de rotina do Air FCA .



8.3.7 Remover a opção DiTi (FCA)

Para remover a opção DiTi, proceda do seguinte modo:

- ✓ Chave para cones DiTi
- 1. Desligue o instrumento.
- 2. Abra o painel de segurança frontal.
- 3. Levante manualmente todas as rodas Z até à posição mais elevada.
- 4. Desloque todas as rodas Z no sentido frontal do instrumento.
- 5. Afaste o mais possível as rodas Z.
- 6. Segure o adaptador de pontas (D) e o tubo de ejeção de pontas (C).
- 7. Desenrosque o cone DiTi (A) com a chave para cones DiTi (B).



8. Puxe o cone DiTi para baixo com cuidado.

8.3.8 Instalar a opção DiTi (FCA)

Para instalar a opção DiTi, proceda do seguinte modo:

- 1. Enrosque o cilindro adaptador (B) no tubo de ejeção de pontas (C).
- 2. Enrosque o cone DiTi (A) no cilindro adaptador. Utilize a chave para cones DiTi.







3. Empurre a tubagem na agulha de plástico até que a tubagem fique bem fixa à opção DiTi.

- 4. Segure o adaptador de pontas (D) e o tubo de ejeção de pontas (C).
- 5. Enrosque o cone DiTi (A) com a chave para cones DiTi (B).



8.3.9 Verificar as pontas fixas

Para verificar as pontas fixas, proceda do seguinte modo:

AVISO

Imprecisões na pipetagem e erros na deteção do líquido!

As pontas dobradas ou o revestimento da ponta danificado causam imprecisões na pipetagem e erros na deteção do líquido.

• Nunca trabalhe com pontas danificadas ou dobradas.



- 1. Desligue o instrumento.
- 2. Abra o painel de segurança frontal.
- 3. Inspecionar as pontas fixas.
- 4. Inspecione o revestimento da ponta fixa com um espelho.

Certifique-se de que as pontas fixas não estão dobradas. Se o revestimento da ponta fixa estiver danificado ou a ponta fixa estiver dobrada, substitua a ponta imediatamente. Consulte a secção "Remover as pontas fixas" [▶ 188].

8.3.10 Remover as pontas fixas

Para remover pontas fixas, proceda do seguinte modo:

- ✓ As pontas fixas foram limpas. Consulte a secção "Tabelas de cuidados do sistema" [▶ 130].
- ✓ As pontas fixas foram verificadas. Consulte a secção "Verificar as pontas fixas" [▶ 187].



- E Tubagem da pipetagem
- 1. Desligue o instrumento.
- 2. Abra o painel de segurança frontal.
- 3. Levante manualmente todas as rodas Z (A) até à posição mais elevada.
- 4. Afaste o mais possível as rodas Z.
- 5. Se estiver instalada uma ponta fixa ajustável, desaperte os quatro parafusos de ajuste da ponta.
- 6. Desenrosque a porca de fixação (C) segurando imediatamente com a outra mão a ponta fixa por baixo da porca de fixação.
- Remova a porca de fixação (C), deslizando-a ao longo do eixo da ponta.
 Evite o contacto entre a porca de fixação e o revestimento da ponta.



- 8. Se a ponta (D) for ajustável, vire a porca de fixação (C) para baixo numa superfície limpa e retire o anel em O e a anilha.
- Se o canal estiver equipado com a opção de volume baixo, desenrosque o rebordo na parte superior da válvula magnética para libertar a tubagem de pipetagem (E) que passa pela roda Z (A).
- 10. Extraia a tubagem de pipetagem (E) uma determinada distância (a) para fora do adaptador de pontas (B) puxando a ponta (D).

Utilize uma lixa de esmeril seca para uma maior aderência na tubagem de pipetagem – não na ponta.

8.3.11 Instalar pontas fixas



Para instalar pontas fixas, proceda do seguinte modo:

1. Puxe cuidadosamente a tubagem de pipetagem aprox. 25 mm (1 pol.) (a) para fora do adaptador de pontas.

Utilize um pedaço de lixa de esmeril para agarrar a tubagem perto da extremidade para assegurar uma maior aderência.

Se, anteriormente, tiver sido instalada uma ponta, corte aprox. 5 mm (0,2 pol.) (b) da tubagem de pipetagem, utilizando uma faca afiada para obter um corte direito.





- **F** Lixa de esmeril
- H Anilha, branca (FEP)
- J Porca de fixação ajustável
- G Alargador de tubagem Te-PS
 - I Anel em O, preto
 - K Parafuso de ajuste da ponta
- 2. Em caso de pontas Te-PS ou pontas de volume baixo:

Utilize o alargador de tubagem Te-PS (G) para alargar a extremidade da tubagem empurrando o alargador de tubagem Te-PS até ao punho para dentro da tubulação enquanto roda a ferramenta.

Enquanto a tubagem ainda está alargada, empurre a ponta Te-PS para dentro extremidade da tubagem em aprox. 4 mm (0,16 pol.).

3. Introduza a porca de fixação na ponta.

Se a ponta for ajustável (por ex., Te-PS), deslize a porca de fixação sobre a anilha (H) e o anel em O (I).

AVISO! Evite o contacto com a extremidade delicada da ponta e o seu revestimento.

- 4. Introduza a ponta e a tubagem de pipetagem no adaptador de pontas.
- 5. Aparafuse a porca de fixação no adaptador de pontas e aperte.

Se a ponta for ajustável (por ex., Te-PS), aperte a porca de fixação de forma que os quatro parafusos de ajuste da ponta (K) fiquem a um ângulo de 45° em relação ao sistema de coordenadas X/Y da mesa de trabalho.

- 6. Limpe as pontas fixas. Consulte a secção Final do dia.
- 7. Execute um teste de precisão da pipetagem conforme definido pelo operador principal.





8.3.12 Remover pontas de perfuração

Para remover uma ponta de perfuração, proceda do seguinte modo:

- ✓ O instrumento está desligado.
- 1. Abra o painel de segurança frontal.
- 2. Levante manualmente todas as rodas Z até à posição mais elevada.
- 3. Desloque todas as rodas Z no sentido frontal do instrumento.
- 4. Afaste o mais possível as rodas Z.





5. Cubra as pontas de perfuração com proteções de pontas de perfuração. Comece com a ponta de perfuração mais recuada.

6. Desenrosque a porca de fixação segurando imediatamente com a outra mão a ponta de perfuração por baixo da porca de fixação.



7. Puxe a tubagem de pipetagem aprox. 25 mm para fora do adaptador de pontas puxando a ponta. Segure a ponta de perfuração na sua parte superior enquanto puxa.



- 8. Puxe a ponta de perfuração para fora da tubagem segurando a tubagem com a outra mão.
- 9. Não remova a proteção da ponta de perfuração. Descarte-a com a ponta de perfuração num recipiente para resíduos biológicos.







8.3.13 Instalar pontas de perfuração

✓ O instrumento está desligado.

- ✓ Está disponível um operador principal.
- 1. Abra o painel de segurança frontal.
- 2. Levante manualmente todas as rodas Z até à posição mais elevada.
- 3. Desloque todas as rodas Z no sentido frontal do instrumento.
- 4. Afaste o mais possível as rodas Z.





 Abra a embalagem da ponta de perfuração. Não remova a proteção da ponta de perfuração (F).

Ordem de instalação das pontas de perfuração: de trás para a frente

- 6. Puxe cuidadosamente a tubagem de pipetagem aprox. 25 mm para fora do adaptador de pontas.
- 7. Empurre a extremidade lisa e cónica da ponta de perfuração para a extremidade da tubagem.



8. Introduza a ponta de perfuração e a tubagem de pipetagem no adaptador de pontas.





9. Enrosque a porca de fixação no adaptador de pontas e aperte manualmente.

- 10. Abra ligeiramente a porca de fixação. Mova a proteção da ponta ligeiramente para baixo para aceder ao eixo da ponta de perfuração. Não remova logo toda a proteção da ponta.
- 11. Rode a ponta de perfuração até que a abertura da ponta fique virada para a parte frontal do instrumento. Mantenha a ponta de perfuração nesta orientação com uma mão e aperte a porca de fixação com a outra mão.
- 12. Verifique se todas as aberturas das pontas estão viradas para a parte frontal do instrumento.







13. Depois de instalar todas as pontas de perfuração, remova todas as respetivas proteções. Comece com a ponta de perfuração mais recuada.

- 14. Contacte um operador principal para repor o contador no FluentControl.
- 15. Contacte um operador principal para realizar um teste do kit de CQ. Consulte os Documentos de referência.
- 16. Execute o método de Fugas do FCA de perfuração .
- 17. Execute um teste de precisão da pipetagem (recomendação: utilize o kit de CQ), conforme definido pelo operador principal.

1



8.3.14 Recolher pontas de perfuração presas

Na eventualidade de uma ponta de perfuração ficar presa e não ser possível recolhê-la através dos comandos de software, esta tem de ser removida manualmente.

Para recolher pontas de perfuração presas, proceda do seguinte modo:

- ✓ O instrumento está desligado.
- 1. Abra o painel de segurança frontal.
- 2. Levante manualmente todas as rodas Z recolhidas até à posição mais elevada.





3. Cubra todas as pontas de perfuração recolhidas com as proteções de pontas de perfuração (A). Comece com a ponta de perfuração mais recuada.



4. Coloque a ferramenta de remoção de pontas de perfuração próxima da ponta presa numa superfície robusta e estável e posicione-a sob a porca de fixação.





- 5. Rode a maçaneta da ferramenta de remoção de pontas de perfuração até que a ponta seja totalmente recolhida.
- 6. Rode a maçaneta na direção oposta e baixe o retrator cerca de 1 cm.
- 7. Remova a ferramenta de remoção da ponta de perfuração.

A ponta de perfuração presa é agora recolhida.

- 8. Limpe a ferramenta de remoção de pontas de perfuração com álcool.
- 9. Verifique se existem danos na ponta de perfuração (por ex., ponta de perfuração dobrada, ponta danificada).
- Substitua a ponta de perfuração caso esteja danificada. Consulte a secção "Remover pontas de perfuração" [▶ 191] e a secção "Instalar pontas de perfuração" [▶ 194].
- 11. Remova todas as proteções das pontas de perfuração segurando na porca de fixação com uma mão e removendo as proteções com a outra mão. Comece com a ponta de perfuração mais recuada.
- 12. Limpe a ferramenta de remoção de pontas de perfuração com álcool.

8.3.15 Verificar o alinhamento do dedo do manipulador

Pode ser necessário o realinhamento dos dedos do manipulador após uma colisão ou ao implementar dedos do manipulador sobresselentes. Isto aplica-se a todos os braços que utilizam manipuladores equipados com dedos do manipulador.



Alinhamento incorreto após uma colisão:

- Analisar a situação.
- Avaliar potenciais causas da colisão ou a causa do desalinhamento do dedo, como uma gaveta desalinhada de um leitor ou dispositivo de lavagem, hotel ou outro segmento programado/posicionado incorretamente.
- Selecionar um procedimento abaixo com base nos requisitos de precisão.
- Se os dedos do manipulador não precisarem de satisfazer requisitos de precisão acima da média, realizar um alinhamento básico. Consultar a secção "Alinhamento básico dos dedos do manipulador para dedos do manipulador do FES" [> 201] ou a secção "Alinhamento básico dos dedos do manipulador para dedos do manipulador fixos" [> 202].



- Caso os dedos do manipulador precisem de satisfazer requisitos avançados (desvio Z <±0,2 mm), realizar o procedimento de alinhamento avançado dos dedos do manipulador. Consultar a secção "Alinhamento avançado dos dedos do manipulador para dedos do manipulador do FES" [▶ 203] ou "Alinhamento avançado dos dedos do manipulador para dedos do manipulador fixos" [▶ 204].
- 3. Os dedos do manipulador podem ser montados com dois parafusos diferentes: a) parafuso Torx M4x12, apertado com um binário de 3 Nm.



 b) parafuso Allen M4x12 em combinação com um anilha tensora (observar a posição de acordo com a ilustração abaixo), apertado com um binário de 3,5 Nm.





Se não estiver disponível uma chave dinamométrica, apertar o parafuso até a anilha ficar plana e a resistência aumentar. Em seguida, apertar ^{1/12} de volta adicional. Tal corresponde aproximadamente a 3,5 Nm.

8.3.16 Alinhamento básico dos dedos do manipulador para dedos do manipulador do FES

Isto aplica-se a todos os braços que utilizam manipuladores equipados com dedos do manipulador.

Para o alinhamento básico, proceder do seguinte modo:

- ✓ O alinhamento incorreto é claramente visível.
- ✓ Não é necessária uma precisão acima da média.
- Chave dinamométrica disponível (com opção de 3 ou 3,5 Nm). Se não estiver disponível uma chave dinamométrica: Parafuso Torx: apertar os parafusos firmemente, mas sem empregar força excessiva.

Parafuso Allen: consultar "Verificar o alinhamento do dedo do manipulador" [▶ 200].



- 1. Remover o dedo do manipulador da cabeça do manipulador.
- 2. Desapertar o parafuso entre o dedo do manipulador e o adaptador do dedo do FES.
- 3. Pressionar o dedo do manipulador contra os batentes superior e traseiro do adaptador conforme apresentado na ilustração abaixo e apertar o parafuso com uma chave dinamométrica (3 ou 3,5 Nm).



8.3.17 Alinhamento básico dos dedos do manipulador para dedos do manipulador fixos

Para o alinhamento básico, proceda do seguinte modo:

- ✓ O alinhamento incorreto é claramente visível.
- ✓ Não é necessária uma precisão acima da média.
- ✓ Chave de fendas dinamométrica disponível (com opção de 3 ou 3,5 Nm).



1. Desaperte o parafuso de fixação (A).



2. Pressione o dedo do manipulador contra os batentes superior e traseiro da montagem da cabeça do manipulador e aperte o parafuso com uma chave de fendas dinamométrica (3 ou 3,5 Nm).

8.3.18 Alinhamento avançado dos dedos do manipulador para dedos do manipulador do FES

Isto aplica-se a todos os braços que utilizam manipuladores equipados com dedos do manipulador.

- ✓ Chave dinamométrica disponível (com opção de 3 ou 3,5 Nm). Se não estiver disponível uma chave dinamométrica: Parafuso Torx: apertar os parafusos firmemente, mas sem empregar força excessiva. Parafuso Allen: consultar "Verificar o alinhamento do dedo do manipulador" [▶ 200].
- 1. Utilizar a ferramenta de deslocação para mudar a altura Z para cerca de 3 mm acima da mesa de trabalho.





2. Desapertar o parafuso entre o dedo do manipulador e o adaptador do dedo do FES.

3. Garantir que o adaptador do dedo do FES está firmemente conectado à cabeça do manipulador.

Os adaptadores do dedo são mantidos no seu lugar numa direção por um *íman.*

 Utilizar a ferramenta de deslocação para mover a altura Z para 0 mm acima da mesa de trabalho. Avançar lentamente até ao último décimo de um milímetro.

Nota: Em caso de dúvidas sobre como aceder ou operar a ferramenta de deslocação, contactar o operador principal.

- 5. Pressionar o dedo do manipulador contra a cabeça do manipulador e a superfície de referência, conforme apresentado na ilustração, e apertar o parafuso com um binário de 3 ou 3,5 Nm.
- Verificar o ajuste ao rodar manualmente a cabeça 90°, 180°, 270°. Um alinhamento incorreto em diferentes posições é indicativo de um alinhamento incorreto da cabeça ou do braço. Neste caso, um FSE tem de verificar o alinhamento.

8.3.19 Alinhamento avançado dos dedos do manipulador para dedos do manipulador fixos

- ✓ Chave de fendas dinamométrica disponível (com opção de 3 ou 3,5 Nm).
- 1. Utilize a ferramenta de deslocação para deslocar a altura Z para cerca de 3 mm.
- 2. Desaperte o parafuso entre o dedo do manipulador e a cabeça do manipulador.
- Utilize a ferramenta de deslocação para mover a altura Z para 0 mm. Desloque lentamente até ao último décimo de um milímetro. Nota: Se não souber como aceder ou operar a ferramenta de deslocação, contacte o seu operador principal.



- 4. Pressione o dedo do manipulador contra a cabeça do manipulador e a superfície de referência e, em seguida, aperte o parafuso com uma chave de fendas dinamométrica de 3 ou 3,5 Nm.
- 5. Verifique o ajuste ao rodar manualmente a cabeça 90°, 180°, 270°. Um alinhamento incorreto em diferentes posições é indicativo de um alinhamento incorreto da cabeça ou do braço. Neste caso, um TASS precisa de verificar o alinhamento.



8.3.20 Verificar o segmento

Verifique se o segmento está fechado.



Fig. 65: Segmento fechado



Fig. 66: Segmento aberto



8.3.21 Remover os pinos de posicionamento

Para remover os pinos de posicionamento, proceda do seguinte modo:

AVISO

Embate ou erro de processo!

Os embates ou outros erros de processamento podem resultar do posicionamento impreciso dos elementos num segmento da plataforma devido a pinos de posicionamento soltos.

- Não opere o Fluent se os pinos de posicionamento estiverem em falta.
- ✓ Os pinos de posicionamento estão partidos.
- 1. Deslize a ferramenta de remoção de pinos sobre o pino de posicionamento.







2. Levante a pega da ferramenta de remoção de pinos e retire o pino do segmento da mesa de trabalho.

8.3.22 Substituir os pinos de bloqueio e os pinos de posicionamento

AVISO

Embate ou erro de processo!

Os embates ou outros erros de processamento podem resultar do posicionamento impreciso dos elementos num segmento da plataforma devido a pinos de posicionamento soltos.

- Não opere o Fluent se os pinos de posicionamento estiverem em falta.
- Os pinos de posicionamento indicados na tabela de cuidados do sistema estão disponíveis.
- 1. Remova o segmento da mesa de trabalho: *Consulte a secção .*
- 2. Pressione o novo pino de bloqueio (B) no orifício (A).





3. Pressione o novo pino de posicionamento (C) no orifício (A).

4. Coloque o segmento na mesa de trabalho: *Consulte a secção .*



9 Embalar, desempacotar, transportar, armazenar e eliminar

Este capítulo inclui informações reguladoras sobre a reciclagem e os rótulos das embalagens que devem ser seguidas.

AVISO

Previna os danos causados por pessoal não qualificado e não autorizado!

Apenas o pessoal da Tecan ou pessoal autorizado pela mesma deve realizar as operações de embalar, desempacotar, transportar e armazenar!

• Consulte o "Apoio ao cliente" [> 213].

Para obter mais informações sobre o deslocamento do instrumento, consulte a secção "Deslocar o instrumento num armário dentro do laboratório" [▶ 139].

9.1 Rótulos das embalagens

A marcação correta e completa das embalagens ajuda a prevenir o manuseamento incorreto, os acidentes, a entrega incorreta, a perda de peso e os danos durante o armazenamento.

Símbolo	Significado	Descrição
	Reciclar	O material da embalagem pode ser reciclado. Não o elimine com o lixo doméstico. As informações sobre o material utilizado para esta embalagem são disponibilizadas abaixo do símbolo.
<u>11</u>	Este lado para cima	Certifique-se de que a embalagem é transportada e armazenada com a face superior, indicada pelas setas, virada para cima. Não a vire.
	Manter seco	Certifique-se de que a embalagem não se molha durante o transporte e o armazenamento.
	Frágil	Manuseie a embalagem com cuidado. Contém produtos frágeis no interior.
*	Proteger da luz solar	Certifique-se de que a embalagem não será exposta ao calor durante o transporte e o armazenamento. Proteja-a de luz solar forte.
		·

Tab. 51: Símbolos da embalagem



Símbolo	Significado	Descrição
	Não empilhar	Não empilhe embalagens. A embalagem não foi concebida para transportar peso extra.

9.2 Eliminação

Esta secção inclui informações regulamentares sobre reciclagem que devem ser seguidas.

AVISO

Reciclagem de acordo com os regulamentos legais aplicáveis!

Cumpra as leis relativas à reciclagem aplicáveis no seu país.

9.2.1 Requisitos locais da União Europeia

A Comissão Europeia adotou a Diretiva relativa aos resíduos de equipamentos elétricos e eletrónicos (REEE; 2012/19/UE).

Desde agosto de 2005, os fabricantes têm a responsabilidade de recuperar e reciclar o equipamento elétrico e eletrónico.

Marcação	Explicação
	 Impactos ambientais negativos associados ao tratamento dos resíduos. Não trate o equipamento elétrico e eletrónico como resíduos urbanos não triados. Separe os resíduos de equipamentos elétricos e eletrónicos.
	 residuos urbanos nao triados. Separe os resíduos de equipamentos elétricos e eletrónicos.

9.2.2 Requisitos locais da República Popular da China

Marcação para a restrição do uso de substâncias perigosas em equipamentos elétricos e eletrónicos

A norma da indústria eletrónica SJ/T11364-2014 da República Popular da China **Marcação para a restrição do uso de substâncias perigosas em equipamentos elétricos e eletrónicos** exige a marcação para a restrição do uso de substâncias perigosas em equipamentos elétricos e eletrónicos.



De acordo com os requisitos especificados na SJ/T11364-2014, todos os produtos elétricos e eletrónicos da Tecan vendidos na República Popular da China são etiquetados com uma marcação para a restrição do uso de substâncias perigosas.

Marcação	Explicação
25	Esta marcação indica que este produto eletrónico contém determinadas substâncias perigosas e pode ser utilizado com segurança durante o período de utilização sem agressão do ambiente, mas deverá entrar no sistema de reciclagem após este período.

9.2.3 Outros requisitos

Marcação	Explicação
Hg	 Esta lâmpada contém mercúrio Recicle-a ou elimine-a conforme exigido pelas leis locais aplicáveis.



10 Apoio ao cliente

Este capítulo explica que ficheiros e informações a Tecan precisa para realizar uma primeira avaliação de um problema.

10.1 Contactos

Contacte o seu distribuidor ou importador local ou um dos seguintes endereços.

Consulte também a nossa página na Internet: www.tecan.com

Tab. 52: Contactos do apoio ao cliente

País/Região	Endereço	Telefone/Fax/	E-mail
Austrália Nova Zelândia Ilhas do Pacífico	Tecan Australia Pty Ltd Unit 2, 475 Blackburn Road Mount Waverly VIC 3149 Austrália	Telefone Telefone Fax E-mail	Número gratuito: 1300 808 403 +61 3 9647 4100 +61 3 9647 4199 helpdesk-aus@tecan.com
Áustria	Tecan Sales Austria GmbH Untersbergstrasse 1a 5082 Grödig Áustria	Telefone Fax E-mail	+43 6246 8933 256 +43 6246 72770 helpdesk-at@tecan.com
Bélgica	Tecan Benelux B.V.B.A. Mechelen Campus Schaliënhoevedreef 20A 2800 Mechelen Bélgica	Telefone Fax E-mail	+32 15 42 13 19 +32 15 42 16 12 tecan-be@tecan.com
Dinamarca	Tecan Denmark, Filial af Tecan Nordic AB, Sverige Lejrvej 29 3500 Værløse Dinamarca	Telefone E-mail	+46 8 7503940 info-dk@tecan.com
França	Tecan France S.A.S.U Tour Swiss Life 1 bd Marius Vivier Merle F- 69 003 Lyon França	Telefone Fax E-mail	+33 4 72 76 04 80 +33 4 72 76 04 99 helpdesk-fr@tecan.com
Alemanha	Tecan Deutschland GmbH Werner-von-Siemens-Straße 23 74564 Crailsheim Alemanha	Telefone Fax E-mail	+49 1805 8322 633 ou +49 1805 TECAN DE +49 7951 9417 92 helpdesk-de@tecan.com



País/Região	Endereço	Telefone/Fax/E-mail	
Itália	Tecan Italia, S.r.l. Via Brescia, 39 20063 Cernusco Sul Naviglio (MI) Itália	Telefone Fax E-mail	+39 800 11 22 91 +39 (02) 92 72 90 47 helpdesk-it@tecan.com
Países Baixos	Tecan Benelux B.V.B.A. Industrieweg 30 NL-4283 GZ Giessen Países Baixos	Telefone Fax E-mail	+31 20 708 4773 +31 183 44 80 67 helpdesk.benelux @tecan.com
Escandinávia	Tecan Nordic AB Sveavägen 159, 1tr SE-113 46 Stockholm Suécia	Telefone Fax E-mail	+46 8 750 39 40 +46 8 750 39 56 info@tecan.se
Espanha Portugal	Tecan Ibérica Instrumentación S.L. C/ Lepanto 151 Bajos E-08013 Barcelona Espanha	Telefone E-mail	+34 93 595 25 31 helpdesk-sp@tecan.com
Suíça	Tecan Schweiz AG Seestrasse 103 8708 Männedorf Suíça	Telefone Fax E-mail	+41 44 922 82 82 +41 44 922 89 23 helpdesk-ch@tecan.com
Reino Unido	Tecan UK Ltd. Theale Court 11-13 High Street Theale, Reading, RG7 5AH Reino Unido	Telefone Fax E-mail	+44 118 930 0300 +44 118 930 5671 helpdesk-uk@tecan.com



Abreviaturas

ADT

Tecnologia de deslocamento de ar

Air FCA

Flexible Channel Arm com sistema de ar

BPL

Boas práticas laboratoriais

CE

Conformité Européenne

CEI

Comissão Eletrotécnica Internacional

CEM

Compatibilidade eletromagnética

cLLD

Deteção capacitiva do nível do líquido

DiTi

Ponta descartável

ΕN

Norma Europeia (European Norm)

FCA

Flexible Channel Arm

FES

Sistema de troca de dedos

FSE

Técnico de assistência local

HEPA

Retenção de partículas de alta eficácia (High-Efficiency Particulate Arrestance)

ISO

Organização Internacional de Normalização (International Organization for Standardization)

LED

Díodo emissor de luz

Liquid FCA

Flexible Channel Arm com sistema de líquido

MCA

Multiple Channel Arm

MCA

Multiple Channel Arm

MCH

Cabeça com múltiplos canais

MET

Marca registada da Euroffins EE enquanto laboratório de testes reconhecido a nível nacional

MIO

Opção de incubadoras monitorizadas (Monitored Incubators Option)

MP

Microplaca

MSA

Manual de software de aplicação

MU

Manual de utilização

NRTL

Laboratório de testes reconhecido a nível nacional



NS

Número de série

PC

Computador pessoal

PP

Polipropileno

QI

Qualificação da instalação

QO

Qualificação operacional

rcf

força centrífuga relativa

RF

Radiofrequência

RGA

Robotic Gripper Arm

RGA Z longo

Braço robótico da garra (Robotic Gripper Arm) longo

RGA Z standard

Braço robótico da garra (Robotic Gripper Arm) de altura standard

RUO

Utilização apenas em investigação

RWP

Bomba de lavagem rápida (Rapid Wash Pump)

SNC

Sistema de Notificação Comum

Te-Shake

Agitador Tecan

Te-VacS

Separador por filtração a vácuo Tecan

USB

Barramento série universal

WEEE

Resíduos de equipamentos elétricos e eletrónicos (Waste Electrical and Electronic Equipment)