

Operating Manual

Fluent®



Název:	Návod k obsluze Fluent			Číslo dílu:	30199937.03
ID:	399706, cz, V2.7	,		Přeloženo z:	-
Verze:	Revize:	Vydání: Historie dokur		entu:	
1	0	30. 8. 2017	První vydání		
1	1	18. 9. 2017	2.3 Rizika aplikad	ce	
1	2	7. 11. 2017	1.3 Zamýšlené p	oužití	
1	3	5. 3. 2018	1.3 Zamýšlené p	oužití	
1	4	16. 4. 2018	3.3 Rozměry a hi	notnosti	
1	5	20. 7. 2018	4.6.2 Upínadlo F	CA	
1	6	26. 2. 2019	2.2 Obecné bezp	ečnostní informac	e
1	7	22. 10. 2020	Informace o míchání a propichování doplněné do částí 1.8.3, 2.5, 2.8.1, 4.6, 4.6.3, 6.6, 7.4.1, 7.4.2 7.5.4, 8.2.3, 8.3.13–15. Informace o rotátoru zkumavek doplněné do částí 4.6.3, 6.5.3, 7.4.1, 7.4.2, 7.5.5, 7.5.6, 8.2.3, 8.3.14. Byly přidány čás 4.6.4 a 7.5.24. Byly aktualizovány části 2.5, 2.9, 2.10, 4.2.2, 4.6, 6.3.5, 7.1, 7.4, 7.5.16, 8.2.4 a 8.3.13–20.		ní doplněné do 6, 7.4.1, 7.4.2, o rotátoru 6.5.3, 7.4.1, /ly přidány části části 2.5, 2.9, .16, 8.2.4
1	8	1. 12. 2020	Byly aktualizovár	ny části 2.2 a 2.8.	
1	9	15. 2. 2021	Byly aktualizovány části 2.7.2, 7.1, 7.2, 7.4.2, 7.5.15, 7.5.24 a 8.2.3 a přidány části 2.7.5, 4.6.4 a 7.5.25.		7.2, 7.4.2, sti 2.7.5, 4.6.4
1	10	25. 11. 2021	Byly aktualizovár	ny části 3.7, 6.5.1.	
2	0	2. 3. 2022	Byla aktualizovár	na část 3.1.	
2	1	14. 4. 2022	Byly přidány část 6.5.1.	i 5.2.5, 6.6, byla a	ktualizována část
2	2	16. 11. 2022	Byla přidána čás a 3.4.	t 4.6.5, byly aktual	izovány části 2.2
2	3	15. 3. 2023	Byly aktualizovár	ny části 1.8.2, 2.4,	3.4 a 7.4.2.
2	4	30. 3. 2023	Byly aktualizovár 4.3.3, 4.3.4, 4.3.5	ny části 2.4, 2.8.1. 5.3, 7.4.2 a 7.4.4.	1, 3.3, 4.3.2,
2	5	16. 1. 2024	Byly aktualizovár 7.2.1, 7.2.2, 7.4.2	ny části 1.9, 2.2, 2 2, 7.5.14, 8.2.2 a 8	.7.1, 2.8.1, 6.2, 7, 3.3.6.

© 2024, Tecan Trading AG, Switzerland, all rights reserved.

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

2	6	6. 6. 2024	Byly aktualizovány části 1, 1.8.4, 2.2, 2.4, 3.7, 4.6, 6.1, byla přidána část 4.6.6.
2	7	5. 12. 2024	Byly aktualizovány části 1.8.4, 3.7, 4.6, 6.5, 7.4.1.2, 7.5.18, 10.1, byla přidána část 2.12.

© 2024, Tecan Trading AG, Switzerland, all rights reserved.

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

Obsah

1	O tomto návodu 8					
	1.1	Rozsah této příručky	8			
	1.2	Výrobce	8			
	1.3	Zamýšlené použití	8			
	1.4	Oblast použití / oblast aplikace	9			
	1.5	Nesprávné použití	9			
	1.6	Záruka	9			
	1.7	Obchodní známky	9			
	1.8	Referenční dokumenty	9			
	1.9	Splnění zákonů a norem	10			
	1.10	Konvence dokumentu	10			
2	Bezp	ečnost	12			
	2.1	Konvence bezpečnostních sdělení	12			
	2.2	Obecné bezpečnostní informace	14			
	2.3	Prohlášení o ochraně osobních údajů pro kameru	15			
	2.4	Rizika aplikace	16			
	2.5	Provozní společnost	25			
	2.6	Validace metody a procesu	25			
	2.7	Kvalifikace uživatele	26			
	2.8	Bezpečnostní prvky	27			
	2.9	Bezpečnostní značky produktu	33			
	2.10	Laserové záření	38			
	2.11	Prohlášení o dekontaminaci	39			
	2.12	Hlášení incidentu	39			
3	Tech	nické údaje	40			
	3.1	Typový štítek	40			
	3.2	Štítek se sériovým číslem	41			
	3.3	Rozměry a hmotnost	42			



	3.4	Napájení	43
	3.5	Datová a napájecí připojení	44
	3.6	Podmínky prostředí	44
	3.7	Emise a odolnost	45
4	Popis	s funkce	47
	4.1	Přehled	47
	4.2	Plošina	47
	4.3	Robotická ramena	51
	4.4	Kapalinový systém (FCA s kapalinovým systémem)	57
	4.5	Mycí systém (MCA 384)	58
	4.6	Doplňky a zařízení	58
5	Řídic	í prvkv	69
•	5.1	Ovládací prvkv	69
	5.2	Uživatelské rozhraní	70
	5.3	Chybové signály a stav přístroje	75
		, , , ,	
	5.4	Stavové LED kontrolky Fluent ID	77
	5.4	Stavové LED kontrolky Fluent ID	77
6	5.4 Oper	Stavové LED kontrolky Fluent ID	77 78
6	5.4 Oper 6.1	Stavové LED kontrolky Fluent ID ace Bezpečnostní pokyny pro tuto kapitolu	77 78 78
6	5.4Oper6.16.2	Stavové LED kontrolky Fluent ID ace Bezpečnostní pokyny pro tuto kapitolu Provozní režimy	77 78 78 79
6	5.4Oper6.16.26.3	Stavové LED kontrolky Fluent ID	77 78 78 79 80
6	 5.4 Oper 6.1 6.2 6.3 6.4 	Stavové LED kontrolky Fluent ID	77 78 79 80 88
6	 5.4 Oper 6.1 6.2 6.3 6.4 6.5 	Stavové LED kontrolky Fluent ID ace Bezpečnostní pokyny pro tuto kapitolu Provozní režimy Uvedení do provozu Před spuštěním metody Spuštění metody	77 78 78 79 80 88 92
6	 5.4 Oper 6.1 6.2 6.3 6.4 6.5 6.6 	Stavové LED kontrolky Fluent ID ace Bezpečnostní pokyny pro tuto kapitolu Provozní režimy Uvedení do provozu Před spuštěním metody Spuštění metody Provoz DeckCheck	77 78 78 79 80 88 92 104
6	5.4 Oper 6.1 6.2 6.3 6.4 6.5 6.6 6.7	Stavové LED kontrolky Fluent ID ace Bezpečnostní pokyny pro tuto kapitolu Provozní režimy Uvedení do provozu Před spuštěním metody Spuštění metody Provoz DeckCheck	77 78 78 79 80 88 92 104 106
6	5.4 Oper 6.1 6.2 6.3 6.4 6.5 6.6 6.7 6.8	Stavové LED kontrolky Fluent ID	77 78 78 79 80 88 92 104 106 108
6	5.4 Oper 6.1 6.2 6.3 6.4 6.5 6.6 6.7 6.8 Péče	Stavové LED kontrolky Fluent ID	77 78 79 80 88 92 104 106 108 110
6	5.4 Oper 6.1 6.2 6.3 6.4 6.5 6.6 6.7 6.8 Péče 7.1	Stavové LED kontrolky Fluent ID	77 78 79 80 88 92 104 106 108 110
6	5.4 Oper 6.1 6.2 6.3 6.4 6.5 6.6 6.7 6.8 Péče 7.1 7.2	Stavové LED kontrolky Fluent ID	77 78 79 80 88 92 104 106 108 110 1110 1111
6	5.4 Oper 6.1 6.2 6.3 6.4 6.5 6.6 6.7 6.8 Péče 7.1 7.2 7.3	Stavové LED kontrolky Fluent ID	77 78 79 80 88 92 104 106 108 110 1110 1111 112



	7.4	Tabulky péče o systém	114
	7.5	Činnosti péče o systém	123
8	Řeše	ní problémů	152
	8.1	Bezpečnostní pokyny pro tuto kapitolu	152
	8.2	Tabulky pro řešení problémů	152
	8.3	Činnosti při řešení problémů	166
9	Baler	ní, vybalování, přeprava, skladování a likvidace	196
9	Baler 9.1	n í, vybalování, přeprava, skladování a likvidace Etikety na obalu	196 196
9	Baler 9.1 9.2	n í, vybalování, přeprava, skladování a likvidace Etikety na obalu Likvidace	196 196 197
9	Baler 9.1 9.2	ní, vybalování, přeprava, skladování a likvidace Etikety na obalu Likvidace	196 196 197
9 10	Baler 9.1 9.2 Záka	ní, vybalování, přeprava, skladování a likvidace Etikety na obalu Likvidace znická podpora	196 196 197 199
9 10	Baler 9.1 9.2 Záka 10.1	ní, vybalování, přeprava, skladování a likvidace Etikety na obalu Likvidace znická podpora Kontakty	196 196 197 199 199
9 10	Baler 9.1 9.2 Záka 10.1	ní, vybalování, přeprava, skladování a likvidace Etikety na obalu Likvidace znická podpora Kontakty	196 197 197 199 199





1 O tomto návodu

Tento Operating Manual popisuje přístroj Fluent a uvádí všechny informace požadované pro jeho bezpečný provoz a udržování v dobrém pracovním stavu. Tento návod je nutné si pečlivě přečíst předtím, než začnete pracovat s přístrojem Fluent a než jej začnete používat.

Tato kapitola uvádí účel tohoto návodu a specifikuje produkt, který je v něm uváděn. Dále vysvětluje použití symbolů a konvencí a uvádí i další obecné informace.

Tento návod se týká vlastního přístroje Fluent. Důležité informace o submodulech naleznete v konkrétních příručkách k modulům.



Tento Operating Manual neobsahuje žádný popis softwaru. Pokud jde o informace o softwaru, použijte příslušný návod k softwaru. Viz část Referenční dokumenty.

1.1 Rozsah této příručky

Tato příručka je platná pro produkt:

- Fluent 480 (číslo dílu 30042011)
- Fluent 780 (číslo dílu 30042021)
- Fluent 1080 (číslo dílu 30042031)

1.2 Výrobce

Adresa výrobce



Tecan Schweiz AG Seestrasse 103 CH-8708 Männedorf Švýcarsko

1.3 Zamýšlené použití

Tento produkt Fluent je plně automatizovaná platforma pro manipulaci s kapalinami pro obecné laboratorní použití. Je určena pro rutinní laboratorní úkoly, jako je pipetování, manipulace s kapalinami a robotická manipulace s laboratorním zařízením v určitých testovacích postupech.



Pro toto zamýšlené použití je nutný software Fluent Gx Assurance Software.

Některé doplňky od firmy Tecan a zařízení třetích stran, které lze používat se zařízením Fluent, jsou určeny jen pro výzkumné použití (RUO).

Jestliže je v zařízení Fluent integrován doplněk nebo zařízení určené jen pro výzkumné použití, nebo pokud není nainstalovaný Fluent Gx Assurance Software, zamýšlené použití se mění takto:

Fluent je plně automatizovaná platforma pro manipulaci s kapalinami pro výzkumné a průmyslové použití. Je určena pro rutinní laboratorní úkoly, jako je pipetování, manipulace s kapalinami a robotická manipulace s laboratorním



zařízením v určitých testovacích postupech (např. buněčné rozbory, biochemické rozbory a správa sloučenin). Tento přístroj není určen pro použití v klinických diagnostických postupech.

1.4 Oblast použití / oblast aplikace

Fluent je možné používat v řadě laboratorních prostředí v souladu se zamýšleným použitím.

V každém prostředí je každá jednotlivá laboratoř odpovědná za validaci přístroje Fluent společně s konkrétními kapalinami a laboratorním vybavením v rámci pracovních postupů nebo metod laboratoře.

1.5 Nesprávné použití

Nesprávné použití může nepříznivě ovlivnit bezpečnostní koncepci přístroje Fluent.

- Fluent se nesmí používat s doplňky nebo komponentami, které nejsou schválené společností Tecan.
- Přístroj Fluent není odolný proti explozi a neměl by být instalován v místech, kde je nebezpečí exploze.
- Fluent by se neměl používat bez přítomnosti funkčních bezpečnostních zařízení.

1.6 Záruka

Fluent se nesmí používat s komponentami, které nejsou schválené společností Tecan.

Použití neschválených komponent může porušit bezpečnostní koncepci přístroje Fluent.

Použití neschválených komponent by zneplatnilo záruku bezpečnosti a shodu s národními a mezinárodními normami, jak je vyžadováno pro certifikaci NRTL a směrnicemi ES atp.

1.7 Obchodní známky

Názvy produktů, ať již se jedná o registrované nebo neregistrované obchodní známky, zmíněné v tomto návodu jsou reprodukovány výhradně pro účely identifikace a zůstávají exkluzivním vlastnictvím svých příslušných vlastníků. Pro účely zjednodušení nejsou symboly obchodních známek, například [®] a ™, v tomto návodu opakovány.

1.8 Referenční dokumenty

Tato část uvádí seznam dokumentů, které jsou potřebné nebo mohou být užitečné při používání přístroje Fluent.

ID dokumentů uvedená níže jsou jejich hlavní čísla. Neobsahují proto informace o jazyku, verzi dokumentu, ani médiu dokumentu (datové úložné médium, výtisk, soubor ke stažení atd.).



Na základě konfigurace vaší objednávky jsou platné také návody k obsluze doplňkových zařízení.



Zkontrolujte rozsah odpovídajícího dokumentu pro zajištění, že máte správnou verzi.

ID dokumentu neodkazuje na objednací informace. Při zadávání objednávek uveďte číslo na vazbě, na pouzdru CD atd.

1.8.1 Příručky k přístroji

- Návod k obsluze Fluent[®] (ID dokumentu 399706)
- Referenční příručka Fluent[®] (ID dokumentu 399937)

1.8.2 Příručky k softwaru

- Příručka k doplňkovému softwaru Tecan pro sledování vzorků (ID dokumentu 393933)
- Příručka k softwaru FluentControl Application Software (ID dokumentu 399935)
- Příručka k aplikačnímu softwaru Introspect (ID dokumentu 400733)
- Příručka k aplikačnímu softwaru MissionControl (ID dokumentu 401940)

1.8.3 Příručky sady QC

- Aplikační příručka sady QC (ID dokumentu 397069)
- Příručka aplikačního softwaru sady QC (ID dokumentu 397070)

1.8.4 Ostatní referenční dokumenty

- Návod k obsluze karuselu Fluent[®] (ID dokumentu 398350)
- HEPA digestoř (ID dokumentu Caron 70072)
- Příručka k aplikaci Frida Reader™ (ID dokumentu 401882)
- Návod k obsluze Te-Shake™ (ID dokumentu 391496)
- Návod k obsluze Te-VacS™ (ID dokumentu 391236)
- Návod k obsluze zásobníku Fluent[®] (ID dokumentu 398658)
- Návod k obsluze MIO2 (ID dokumentu 394934)
- Návod k obsluze Resolvex i300 (ID dokumentu 402756)

1.9 Splnění zákonů a norem

Na přístroj Fluent se vztahují následující prohlášení a certifikace:

- Prohlášení o shodě ES s platnými směrnicemi EU (značka CE)
- Certifikace celostátně uznávané zkušební laboratoře (NRTL)
- Certifikace schématu CB (IECEE) (značka CB)

Podrobnější informace o značení uvádí část "Typový štítek" [> 40].

1.10 Konvence dokumentu

Křížové odkazy se objevují následujícím způsobem – např.:
Viz část "Bezpečnost" [▶ 12]
"Bezpečnost" odkazuje na odpovídající záhlaví části
Číslo strany je uvedeno v hranatých závorkách

Předpoklady Předpoklady se objevují následujícím způsobem – např.:



✓ Byly přečteny "Obecné bezpečnostní informace".

Тіру

Doplňující rady se objevují následujícím způsobem – např.:



Ohledně bezpečnostních konvencí a symbolů viz kapitolu "Bezpečnost" [> 12].

llustrace

llustrace mohou znázorňovat verze komponent, které nejsou relevantní pro váš produkt Fluent.



2 Bezpečnost

Tato kapitola popisuje bezpečnostní koncepci produktu Fluent, uvádí obecná pravidla správného chování a varování před riziky spojenými s používáním produktu Fluent.

2.1 Konvence bezpečnostních sdělení

2.1.1 Signální slova

Tab. 1: Signální slova

Signální slovo	Význam
A NEBEZPEČÍ	Indikuje nebezpečnou situaci, která, pokud jí nebude zabráněno, povede k úmrtí nebo závažnému poranění.
A VAROVÁNÍ	Indikuje nebezpečnou situaci, která, pokud jí nebude zabráněno, by mohla vést k úmrtí nebo závažnému poranění.
A UPOZORNĚNÍ	Indikuje nebezpečnou situaci, která, pokud jí nebude zabráněno, by mohla vést k malému nebo střednímu poranění.
OZNÁMENÍ	Indikuje situaci, která nesouvisí s nebezpečím, ale pokud jí nebude zabráněno, mohla by vést k poškození nebo nefunkčnosti zařízení, nebo k nesprávným výsledkům procesu.



2.1.2 Bezpečnostní symboly





2.2 Obecné bezpečnostní informace

A VAROVÁNÍ

Přístroj Fluent je zkonstruován a sestaven v souladu se současnou moderní technologií a uznávanými technickými bezpečnostními předpisy. Přesto se mohou vyskytnout rizika ohrožující uživatele, majetek a prostředí, pokud by se přístroj Fluent používal bez náležité péče a pozornosti.

Bezpečnost všech uživatelů a personálu závisí na přísném dodržování těchto bezpečnostních pokynů a na znalosti bezpečnostních varování uvedených v tomto návodu.

- Věnujte prosím velkou pozornost následujícím obecným bezpečnostním informacím.
- Tento návod musí být vždy k dispozici všem osobám, které provádějí úkoly v něm popsané.
- Vždy používejte napájecí kabel dodaný s přístrojem.
- Nepoužívejte napájecí kabel s jinými produkty.
- Je třeba důsledně dodržovat právní předpisy, jako jsou místní, státní a federální zákony týkající se používání a aplikace, taktéž manipulace s nebezpečnými materiály v souvislosti s přístrojem Fluent.
- Provozní společnost je odpovědná za definování pokynů v souladu s firemními postupy a místními právními požadavky. Je potřeba důsledně dodržovat pokyny provozovatele.
- Dodržujte správné podmínky prostředí pro skladování a provoz.
- Konstrukční změny bezpečnostních zařízení jsou zakázány.
- Poškozené bezpečnostní zařízení musí být neprodleně vyměněno, jak je popsáno v tomto návodu.
- Přístroj Fluent nesmí být nijak modifikován bez předchozí konzultace a písemného souhlasu společnosti Tecan. Autorizované modifikace systému může provádět pouze FSE s certifikací pro opravy a inovaci přístroje Fluent. Společnost Tecan odmítne jakoukoli reklamaci vyplývající z neoprávněných modifikací.
- Nebezpečí požáru způsobené nesprávným používáním přístroje Fluent. Přístroj Fluent nesmí být instalován v místech, kde je nebezpečí exploze.
- Nebezpečí požáru způsobené hořlavými kapalinami nebo kapalinou v systému.
- Zabraňte vzniku a hromadění hořlavých výparů.
- S použitými látkami nebo vzorky a činidly zpracovávanými pomocí přístroje Fluent může být spojeno chemické, biologické a radioaktivní nebezpečí (např. při nakládání a vykládání). Totéž platí pro likvidaci odpadu.
 - Vždy mějte na paměti možná nebezpečí spojená s těmito látkami.
 - Používejte vhodný ochranný oděv, ochranné brýle, respirátory a rukavice.
 - Nakládání s látkami a likvidace odpadů mohou podléhat místním, státním nebo federálním zákonům nebo předpisům, které se týkají zdraví, životního prostředí nebo bezpečnosti. Přísně dodržujte příslušná ustanovení.
- Jakékoli znečištění musí být okamžitě odstraněno způsobem popsaným v tomto návodu.



- Uživatel je odpovědný za to, že přístroj Fluent je vždy provozován za správných podmínek a že údržba, servis a opravy jsou prováděny pečlivě, podle plánu a pouze autorizovaným personálem.
- Riziko nesprávných výsledků měření. Po provedené údržbě nebo ošetření systému smí být provoz obnoven až po ověření správných provozních podmínek systému.
- Pro zajištění dobrého výkonu a spolehlivosti systému vždy používejte doporučený spotřební materiál v rámci doby použitelnosti a originální náhradní díly pro údržbu a opravy.
- Při kontaktu pokožky se systémovou kapalinou přístroje by mohlo dojít ke zranění.

- Vždy noste ochranný oděv v souladu s GLP.

- Těžké břemeno! Nezvedejte přístroj.
- Neprovozujte systém bez podnosů a segmentů pracovní plochy.
- Podnosy pracovní plochy zachycují rozlité kapaliny, které se mohou rozlít v prostoru ručního vkládání na pracovní plochu. Systém by měl být provozován s co největším počtem podnosů pracovní plochy nainstalovaných pod pracovní plochou, aby byly zachyceny všechny rozlité kapaliny. Neprovozujte systém bez podnosů pracovní plochy.
- Pokud není povoleno přenášení, důrazně se doporučuje používat jednorázové špičky s filtry.
- Možné rozdrcení. Nepokládejte na pracovní plochu zařízení bez modelových dat Tecan.
- Nástavec 300 je určen pro maximální zatížení 40 kg (88 lb) a pouze pro použití s doplňky, které jsou lehčí než 40 kg (88 lb).
- Produkt Fluent je dodáván s bezpečnostní značkou biologického nebezpečí, kterou by měl uživatel použít v případě použití biologicky nebezpečných látek. Štítek nalepte na přední dveře na místo, které je viditelné uživateli a vhodné k použití. Viz část Bezpečnostní značky produktu.
- Doplňky používané na pracovní desce přístroje Fluent mohou generovat silné magnetické pole, které může narušovat funkci zdravotnických prostředků (ať už implantovaných nebo nošených), jako je kardiostimulátor nebo inzulínová pumpa. Přístroj Fluent je dodává s bezpečnostním štítkem upozorňujícím na silná magnetická pole, který je nutno nalepit na přední dveře v poloze viditelné pro uživatele v případě práce s doplňky generujícími silná magnetická pole.
- Instalaci ethernetového kabelu kamer DeckCheck provede FSE, přičemž kabel musí být za všech okolností připojen k počítači přístroje Fluent (EMC). Rozhraní Ethernet nesmí být připojeno k síti.
- Jen pro obyvatele Kalifornie: Tento výrobek vás může vystavit působení chemických látek, například olovu, které je v Kalifornii považováno za látku, která způsobuje rakovinu, vrozené vady nebo jiné reprodukční poškození. Další informace uvádí www.P65Warnings.ca.gov/product.

2.3 Prohlášení o ochraně osobních údajů pro kameru

Systém Fluent je vybaven kamerami umístěnými na vnitřním předním profilu. Kamery jsou namířeny na plošinu a na zadní plošinu. Přes boční panely z akrylového skla je možný pohled směrem dolů.



- Uživatel je odpovědný za upozornění osob v místnosti, že jsou kamery v provozu.
- Uživatel je odpovědný za to, že na pořízených snímcích nebude možné identifikovat personál, například pokud přístroj (z boku) sousedí s prostorem stolu nebo pokud jsou výřezy v zadním či bočním panelu nebo pokud je místo zadní stěny použit panel z akrylového skla.

2.4 Rizika aplikace

Systémové funkce / modul	Možný režim selhání	Potenciální účinek selhání	Možná/potenciální příčina	Značení nebo zmírnění
Systém	Nedostatečn á údržba	Bezpečnost nebo zdraví uživatelů: Potenciální kontaminace přístroje	Chyba použití: Nedodržení pokynů uvedených v návodu k obsluze nebo pokynů pro údržbu	Uživatel musí zajistit používání vhodného spotřebního materiálu a dodržování pokynů pro preventivní údržbu (viz část Péče o systém). Uživatel musí používat ochranný oděv, rukavice a ochranné brýle v souladu se zásadami GLP a platnými místními předpisy.
Systém	Požár	Bezpečnost nebo zdraví uživatelů: Požár v laboratoři operátorů (hořící přístroje)	Plyn z těkavých hořlavých kapalin; šíření jiskry z elektronické desky	Přístroj není chráněný proti výbuchu a zákazník musí zajistit, aby nedocházelo k vytváření vysoké koncentrace výparů (viz část Obecné bezpečnostní informace).
Modul FCA a Air FCA	Opotřebení mechaniky osy Z (nadprůměrn é opotřebení)	Bezpečnost nebo klinické podmínky vzorku: Potenciálně nesprávné polohování Z v laboratorním vybavení	Nadprůměrné použití přístroje v kombinaci s použitím jednorázových špiček Vysoké procento kroků propichování v aplikaci	Systém informuje uživatele, pokud osa Z dosáhne 90 % očekávané životnosti.
Modul FCA a Air FCA	Opotřebení mechaniky osy P (nadprůměrn é opotřebení)	Bezpečnost nebo klinické podmínky vzorku: Potenciálně nesprávné polohování ve směru osy P v laboratorním vybavení	Nadprůměrné použití přístroje v kombinaci s použitím jednorázových špiček Vysoké procento kroků propichování v aplikaci	Systém informuje uživatele, pokud osa P dosáhne 90 % očekávané životnosti.



Systémové funkce / modul	Možný režim selhání	Potenciální účinek selhání	Možná/potenciální příčina	Značení nebo zmírnění
Modul FCA a Air FCA	Abraze ozubeného kola pohonu X (nadprůměrn é použití)	Bezpečnost nebo klinické podmínky vzorku: Potenciální kontaminace vzorků polyamidovými částicemi	Nadprůměrné použití přístroje v kombinaci s umístěním kritického laboratorního vybavení na zadní straně přístroje	Neumisťujte prvky citlivé na částice (např. vzorky a činidla) na zadní stranu přístroje nebo na horní část laboratorního vybavení umístěte ochranu proti částicím (např. víka).
Modul FCA a Air FCA	Rušivé signály způsobené propíchnutím přepážky	Bezpečnost nebo klinické podmínky vzorku pacienta: nesprávná položka cLLD vedoucí k aspiraci vzduchu a potenciálně falešným výsledkům	Interakce špičky s přepážkou/fólií	Pracujte pouze s nevodivými fóliemi pro propichovací aplikace ve spojení s detekcí hladiny kapaliny na modulech FCA a Air FCA. Viz referenční příručku. Uživatel musí validovat detekci kapaliny v kombinaci s propíchnutím pro moduly FCA a Air FCA.
Modul FCA a Air FCA	Chybná úprava vzorku, chybné cLLD v důsledku pěny nebo bublinek v lahvičce s činidlem	Bezpečnost procesu: nesprávně zpracované vzorky	Bublinky nebo pěna v lahvičce s činidlem způsobují nesprávné cLLD a možnou aspiraci vzduchu modulem FCA nebo Air FCA	Uživatel je odpovědný za validaci aplikace/procesu s ohledem na vhodnou přípravu vzorku.
Modul FCA a Air FCA	Ucpání špičky	Bezpečnost nebo klinické podmínky vzorku pacienta: Potenciálně chybný pipetovaný objem	Aspirace na dně nádoby (zablokování špičky)	Uživatel musí validovat aplikaci, aby zabránil aspiraci příliš blízko úrovně Z-max vlastního laboratorního vybavení.



Systémové funkce / modul	Možný režim selhání	Potenciální účinek selhání	Možná/potenciální příčina	Značení nebo zmírnění
Modul FCA a Air FCA	Hadicový systém FCA: Růst mikroorganis mů	Bezpečnost nebo klinický stav vzorku pacienta: nesprávně pipetovaný objem nebo kontaminace vzorků	Růst mikroorganismů (biologická vrstva na vnitřním povrchu)	Jako systémovou kapalinu pro modul FCA používejte deionizovanou vodu a provádějte denní údržbu zahrnující proplachování systému podle pokynů pro každodenní péči o systém (viz část Péče o systém), přičemž používejte pouze čisticí prostředky povolené pro tento úkon.
Modul MCA 96 nebo MCA 384/96	Přetečení kapaliny vzorku v mikrodesce při pipetování	Bezpečnost nebo klinická podmínka vzorku pacienta: Potenciální křížová kontaminace vzorků (přetečení)	Chybně definované úrovně Z uživatelem (např. aspirace z polohy Z-max)	Definujte bezpečné polohy pro aspiraci a dávkování. Viz referenční příručku.
Modul MCA 96 nebo MCA 384/96	Vzorky zcela nebo částečně minou zamýšlenou polohu v režimu volného dávkování	Bezpečnost nebo klinická podmínka vzorku pacienta: Potenciální křížová kontaminace	Elektrostatické náboje na konci špičky způsobené používáním přístroje mimo stanovené podmínky vedou k tomu, že vzorek zůstává viset na špičce nebo je nekontrolovaně rozstříknut	Uživatel musí dodržovat stanovené provozní podmínky pro manipulaci s kapalinami pro modul MCA, zejména pokyny týkající se minimální požadované vlhkosti (viz část "Podmínky prostředí" [▶ 44]). Uživatel musí nastavit výšku dávkování vždy v jamce. Viz referenční příručku.
Modul MCA 96 nebo MCA 384/96	Míchání vzduchu místo kapaliny (vzorek/ činidlo) pro pipetování směsi	Bezpečnost nebo klinická podmínka vzorku pacienta: Vzorky mohou být nesprávně zpracovány s následkem falešných výsledků	Nevhodné parametry sledování v důsledku nesprávné kombinace špiček a mikrodesek	Uživatel by měl porovnat skutečný a virtuální pracovní stůl pomocí názvu laboratorního vybavení na virtuálním pracovním stole. Uživatel by měl dodržovat jedinečné barevné provedení (specifické pro typ špičky) a značení (pro filtr a bez filtru) boxů DiTi. Uživatel musí před zahájením procesu zkontrolovat rozvržení pracovního stolu.



Systémové funkce / modul	Možný režim selhání	Potenciální účinek selhání	Možná/potenciální příčina	Značení nebo zmírnění
Specifická jednorázová špička	Získání DiTi: Namontovan ý nesprávný typ špičky	Bezpečnost nebo klinická podmínka vzorku pacienta: Potenciálně žádný nebo krátký aspirovaný vzorek Potenciální křížová kontaminace vzorků	Chyba použití: Nesprávné rozložení pracovní plochy: uživatel pokládá box špiček na nesprávné místo: špičky jsou kratší, než se očekává Nesprávné rozložení pracovní plochy: uživatel pokládá box špiček na pracovní stůl s nefiltrovanými špičkami namísto filtrovaných špiček Nesprávné rozložení pracovní plochy: uživatel pokládá box špiček Nesprávné rozložení pracovní plochy: uživatel pokládá box špiček na nesprávné místo: špička má menší objem, než se očekává (např. 100 µl místo 200 µl); délka špičky odpovídá očekávání; kapalina aspirovaná v MCH	Uživatel by měl porovnat skutečný a virtuální pracovní stůl pomocí názvu laboratorního vybavení na virtuálním pracovním stole. Uživatel by měl dodržovat jedinečné barevné provedení (specifické pro typ špičky) a značení (pro filtr a bez filtru) boxů DiTi. Uživatel musí před zahájením procesu zkontrolovat rozvržení pracovního stolu. Mechanická konstrukce zajišťuje viditelnost bílého filtru. Referenční příručka obsahuje informace o barevném značení boxů DiTi, o rozdílu délek a filtrovaných DiTi. Viz referenční příručku.
Specifická jednorázová špička	Neúplné pouštění špiček: Některé kontaminova né špičky zůstávají viset na hlavici a padají na destičky se vzorky	Bezpečnost nebo klinická podmínka vzorku pacienta: Potenciální křížová kontaminace	Způsobené elektrostatické náboje	Uživatel musí dodržovat stanovené provozní podmínky pro manipulaci s kapalinami pro modul MCA, zejména pokyny týkající se minimální požadované vlhkosti (viz část "Podmínky prostředí" [▶ 44]). Jednorázové špičky nejsou určeny k opakovanému použití.



Systémové funkce / modul	Možný režim selhání	Potenciální účinek selhání	Možná/potenciální příčina	Značení nebo zmírnění
Modul RGA	Ztráta desky v důsledku srážky s nesprávně vyrovnaným laboratorním vybavením	Bezpečnost procesu: Ztráta desky, ztráta vzorků	Pokud jsou na sebe naskládány více než 4 mikrodesky, během přepravy může dojít ke špatnému vyrovnání	Pohyby desek musí být před spuštěním skriptů se skutečnými vzorky validovány. Viz kontrolní seznam validace v příručce k aplikačnímu softwaru.
Module FluentContro I Software	WorktableBa se: byl vykázán chybný stav DiTi	Bezpečnost procesu: Křížová kontaminace / chybné výsledky	Křížová kontaminace z důvodu nesprávných informací o stavu použití špiček	Nepoužívejte možnost "Set Tips Back" (Nastavit špičky zpět), pokud režim selhání vede k riziku vysoké závažnosti.
Module FluentContro I Software	Hodnota Core.Scriptin g.Programmi ng SetVariable v runtime: chybná hodnota	Bezpečnost procesu: chybné výsledky	Chyba v softwaru: proměnná je nastavena na chybnou hodnotu	Proveďte validaci aplikace pro specifický zdroj, cíl a rozsahy proměnné. Viz kontrolní seznam validace v příručce k aplikačnímu softwaru.
Module FluentContro I Software	Hodnota Core.Scriptin g.Programmi ng QueryVariabl e v runtime nebo spuštění skriptu: chybná prezentace UI / přijetí hodnoty UI	Bezpečnost procesu: chybné výsledky	Numerická hodnota je v UI nesprávně naformátovaná nebo převedená	Proveďte validaci aplikace pro specifický zdroj, cíl a rozsahy proměnné. Viz kontrolní seznam validace v příručce k aplikačnímu softwaru.
Module FluentContro I Software	Hodnota Core.Scriptin g.Programmi ng ImportVariabl e v runtime: byla importována chybná hodnota	Bezpečnost procesu: chybné výsledky	Ze zdroje importu je vyvolána nesprávná hodnota	Proveďte validaci aplikace pro specifický zdroj, cíl a rozsahy proměnné. Viz kontrolní seznam validace v příručce k aplikačnímu softwaru.



Systémové funkce / modul	Možný režim selhání	Potenciální účinek selhání	Možná/potenciální příčina	Značení nebo zmírnění
Module FluentContro I Software	Hodnota Core.Scriptin g.Programmi ng ExportVariab le v runtime: byla exportována chybná hodnota do souboru	Bezpečnost procesu: chybné výsledky	Do souboru exportu byla zapsána chybná hodnota	Proveďte validaci aplikace pro specifický zdroj, cíl a rozsahy proměnné. Viz kontrolní seznam validace v příručce k aplikačnímu softwaru.
Module FluentContro I Software	API: Získání/ nastavení proměnné nebo vyřešení výrazu selhalo	Bezpečnost procesu: chybné výsledky	Byla vyvolána chybná proměnná nebo přiřazena chybná hodnota / byl vrácen chybný výraz	Proveďte validaci aplikace pro specifický zdroj, cíl a rozsahy proměnné. Viz kontrolní seznam validace v příručce k aplikačnímu softwaru.
Světlo UVC	Nesprávné použití v aplikaci	Nedostatečná účinnost	Nesprávné použití v aplikaci	Viz specifické pokyny v části "Optické záření (UVC)" [▶ 32].
Rotátor zkumavek / propichovací špičky (Míchání a propichová ní)	Nesprávné použití v aplikaci	Nedostatečná účinnost	Nesprávné použití v aplikaci	Viz specifické pokyny v části "Míchání a propichování" [▶ 62].
Čtečka Frida	Nesprávné použití v aplikaci	Nedostatečná účinnost	Nesprávné použití v aplikaci	Viz specifické pokyny v části "Čtečka Frida" [▶ 65].
Jakýkoliv	Neefektivní použití v aplikaci	Nedostatečná účinnost v aplikaci	Nedostatečná péče o systém	Viz pokyny pro jednotlivé díly v kapitole Péče o systém.
Zpracování potenciálně nebezpečnýc h materiálů	Kontaminace potenciálně nebezpečný mi materiály	Potenciální rizika pro uživatele, majetek a životní prostředí.	Nedodržování obecných bezpečnostních informací	Viz pokyny pro jednotlivé díly v části Obecné bezpečnostní informace.



Systémové funkce / modul	Možný režim selhání	Potenciální účinek selhání	Možná/potenciální příčina	Značení nebo zmírnění
MCA 96, cLLD	Nesprávný výsledek měření	Chybné měření: Zjištěná hladina kapaliny se nevztahuje na všechny jamky mikrodesky: Chybný výsledek testu nebo ztráta vzorku	Použití nevhodného/ nekompatibilního laboratorního vybavení	Funkci cLLD lze poté použít pouze ve vaničkách.
MCA 96, obecně	Nekompatibil ita spotřebního materiálu / komponent / modulů	Ztráta DiTi během procesu: Potenciální nenapravitelná ztráta vzorku. Potenciální křížová kontaminace.	Použití nevhodného/ nekompatibilního spotřebního materiálu. Je vydán příkaz k odběru s posuvem, který není kompatibilní s boxem nebo zásobníkem, z kterého mají být instalovány DiTi (např. nesprávná verze zásobníku DiTi). DiTi jsou nesprávně instalovány a během procesu dojde k jejich ztrátě	Použijte boxy DiTi s kombinovaným zásobníkem Tecan.
MCA 96, cLLD	Nesprávný výsledek měření	Chybné měření: Falešně pozitivní detekce subsystému cLLD.	Uživatel umístí na pracovní stůl jiná DiTi, než je uvedeno v softwaru. Počet nasazených DiTi se liší od počtu DiTi očekávaného softwarem z důvodu zásahu uživatele (např. uživatel vyjme některá DiTi). Falešně pozitivní detekce z důvodu nesprávné prahové hodnoty cLLD.	Je důležité, aby počet DiTi použitý pro detekci cLLD odpovídal počtu, který je uveden v softwaru.



Systémové funkce / modul	Možný režim selhání	Potenciální účinek selhání	Možná/potenciální příčina	Značení nebo zmírnění
MCA 96, obecně	Kontaminace vzorku	Kontaminace vzorku v důsledku rozlití po nárazu. Chybný výsledek testu nebo ztráta vzorku.	Nesprávně nastavené parametry (nesprávný vektor pro odběr nebo snížení, nesprávný typ nástroje). Kolize hlavy modulu MCA 96 během vektorového pohybu. Například s vysokým laboratorním vybavením na sousedních soupravách.	Existují rizika spojená s chybným použitím funkce vektorového pohybu.
MCA 96, obecně	Kontaminace vzorku	Bezpečnost nebo klinická podmínka vzorku pacienta: Kontaminace vzorku: Potenciální křížová kontaminace v důsledku oděru pásu upínadla.	Oděr materiálu pásu v průběhu životnosti: Odřené částečky pásu osy G mohou spadnout do vzorků pacientů na pracovním stole a posléze vzorky chemicky kontaminovat.	Odřené částečky (z řemenů) a prach mohou spadnout na pracovní stůl a kontaminovat vzorky nebo chemické látky. Tomu lze zabránit zakrytím laboratorního vybavení obsahujícího citlivé kapaliny umístěním víček.
MCA 96, obecně	Kontaminace vzorku	Bezpečnost nebo klinická podmínka vzorku pacienta: Potenciálně chybné výsledky: Potenciální křížová kontaminace v důsledku rozlití do sousedních jamek na desce.	Rozstříknutí nebo rozlití kapaliny ze vzorku. Kontaminace kuželů DiTi kapalinou obsaženou v uchopeném laboratorním vybavení, například když je plné laboratorní vybavení vyhozeno do odpadu a kapalina se vylije.	Laboratorní vybavení před likvidací nejprve vyprázdněte.



Systémové funkce / modul	Možný režim selhání	Potenciální účinek selhání	Možná/potenciální příčina	Značení nebo zmírnění
MCA 96, obecně	Přenos vzorku nebo činidla (FC)	Potenciální křížová kontaminace vzorků způsobená kontaminovaný mi kužely DiTi a blokem válce.	Zařízení se pokouší zvednout DiTi. Software nedokáže detekovat chybějící DiTi, a proto pokračuje bez signalizace chyby. Kužely DiTi se potenciálně dostanou do kontaktu s kapalinou z vaničky a nasají ji.	Modul MCA 96 nedokáže rozpoznat prázdné boxy pro špičky, pokud jsou používány boxy, u nichž je okraj špičky v rovině s horním povrchem boxu.
MCA 96, obecně	Rozstříknutí nebo rozlití kapaliny ze vzorku	Kontaminace vzorku způsobená rozlitím z důvodu nedostatečně pevného laboratorního vybavení.	Rozstříknutí nebo rozlití kapaliny ze vzorku z důvodu použití nedostatečně pevného laboratorního vybavení.	K přepravě kapalin používejte pouze dostatečně pevné laboratorní vybavení, aby nedocházelo k jeho deformaci.
MCA 96 obecně	Prst upínadla může být poškozen v důsledku nárazu	Poškozený prst upínadla		Zkontrolujte prsty upínadla a v případě poškození je vyměňte.
Mycí stanice, míchání a propichová ní	Kontaminace vzorku	Odpad správně neodtéká. Kontaminace vzorku z důvodu nesprávného vypouštění odpadu a nesprávného mytí špiček.	Nesprávně umyté špičky v důsledku přetečení mycí stanice způsobeného zablokováním spojovacího potrubí k mycí stanici.	Systém vyžaduje pravidelnou údržbu. Přípojky mycí stanice je nutno vždy po 2–3 měsících práce s plnou krví vyměnit, aby nedošlo k jejich zanesení. Vypouštění potenciálně korozivních kapalin, jako je 2% roztok chlorové dezinfekce, skrze mycí stanici, přípojky a hadičky nelze provádět bez následného proplachu neutrální kapalinou (např. vodou), aby nedocházelo ke korozi.



2.5 Provozní společnost

•

Provozovatel musí zajistit, aby přístroj Fluent a zejména bezpečnostní prvky fungovaly správně a aby všichni pracovníci, kteří s přístrojem přicházejí do styku, byli náležitě proškoleni.

Odpovědnosti

- Definování procesů v souladu se standardními provozními postupy.
- Zajištění dokončení instalačních a provozních kvalifikací (IQ OQ).
- Zajištění, aby veškerý personál, který je ve styku s přístrojem Fluent, byl náležitě vyškolen.
- Zajištění dostupnosti vhodného ochranného oděvu a prostředků.
- Zajištění údržby a bezpečného provozu přístroje Fluent.
- Požadavky na dodržování laboratorních bezpečnostních předpisů a směrnic.

2.6 Validace metody a procesu

Validace metody a procesu.

Při provádění validace metod a procesu věnujte pozornost následujícím skutečnostem:

- Pokud používáte fixní špičky s modulem MCA 384 nebo FCA, ujistěte se, že postup mytí je účinný pro očekávaný rozsah koncentrace vzorku a citlivosti testu.
- Zkontrolujte, zda pipetované objemy splňují požadavky na přesnost a správnost automatizovaného procesu.
- Při použití laboratorního vybavení jiného než Tecan nebo vlastního vybavení a při aspiraci se sledováním se ujistěte, že definice nádoby je správná (tj. že je použita vhodná rychlost pro sledování), aby nedošlo k aspiraci vzduchu.
- Funkčnost separátoru fází byla ověřena při použití se standardními jednorázovými špičkami Tecan 1 ml a jednorázovými špičkami Tecan 1 ml se širokým průměrem. Další informace o podporovaném spotřebním materiálu Tecan naleznete v referenční příručce (viz část "Referenční dokumenty" [> 9]).
- Validujte detekci kapaliny na přenosové stanici Fluent Stacker.
- Validujte správné používání mycí stanice MCA v rámci aplikace.
- Validujte aplikaci s ohledem na správné objemy pipetování a sledování.
- Validujte aplikaci za účelem prevence aspirace příliš blízko Z-max vlastního laboratorního vybavení.
- Validujte aplikaci propichování s ohledem na potřebné držáky (aktivní nebo pasivní).
- Pokud chemikálie a laboratorní vybavení nejsou odebrány, je třeba vyhodnotit dopad světla UVC na chemikálie a laboratorní vybavení přítomné na plošině a provést validaci testu.
- Zahrňte manuální kontrolu správných objemů pipetování po spuštění.
- Personál musí být informován o prohlášení o ochraně osobních údajů v souvislosti s kamerami (viz "Prohlášení o ochraně osobních údajů pro kameru" [▶ 15]).

Odpovědnost hlavního operátora



2.7 Kvalifikace uživatele

Pracovníci laboratoře musí být plně kvalifikovaní a vyškolení k obsluze přístroje Fluent. Práce popsané v tomto Operating Manual mohou provádět pouze oprávnění pracovníci s níže uvedenou kvalifikací.

Laboratorní personál musí:

- mít odpovídající technické vzdělání,
- · znát bezpečnostní předpisy a směrnice pro laboratoře,
- znát pokyny k bezpečnostním prvkům přístroje,
- používat ochranný oděv a vybavení,
- znát a dodržovat správnou laboratorní praxi,
- a přečíst si pokyny v návodu k obsluze a rozumět jim.

Společnost Tecan doporučuje, aby operátor navštěvoval školicí kurz operátorů. O dostupných kurzech se informujte v zákaznickém servisu společnosti Tecan. Viz část "Zákaznická podpora" [> 199].

2.7.1 Operátor

Operátor (laborant) pracuje pro provozní společnost.

- Žádné specifické znalosti aplikací nebo systémů
- Znalost místních jazyků
- Znalost angličtiny je výhodou

Operátor má přístupová práva k aplikačnímu softwaru umožňující spouštět metody a provádět péči o systém a získá potřebné školení od hlavního operátora.

2.7.2 Hlavní operátor

Hlavní operátor (specialista aplikace) podporuje provozní společnost nebo pracuje pro stejnou společnost.

Požadované dovednosti

Požadované

dovednosti

- Rozsáhlé znalosti aplikací
- Omezená znalost systému
- Znalost místních jazyků
- Znalost angličtiny
- Důkladná znalost příručky příslušného softwaru

Odpovědnosti

- Pokyny pro operátora
- Psaní, spouštění a validace metod
- Pomoc operátorovi při řešení problémů s přístrojem



2.8 Bezpečnostní prvky

DEPOZORNĚNÍ

Pohyblivé díly

Ochranné a bezpečnostní prvky nainstalované na produktu Fluent nesmí být během provozu odstraněny, vyřazeny z činnosti nebo překonány.

 Pokud jsou některá zařízení demontována (např. kvůli údržbě), musí být před obnovením provozu všechna ochranná a bezpečnostní zařízení znovu nainstalována, aktivována a zkontrolována.

Bezpečnostní panely a bezpečnostní snímače jsou nedílnou součástí produktu Fluent, zatímco zámky dveří přístroje a zámky dveří skříně mohou být součástí pouze některých konfigurací systému.

2.8.1 Bezpečnostní panely

Přístroj Fluent je chráněn bezpečnostními panely:

Přední bezpečnostní panel lze otevřít a je vybaven snímači dveří, které spouštějí aktivní zastavení. Přední bezpečnostní panel lze uzamknout volitelnými zámky dveří.

Zařízení Fluent s rameny MCA 96 nebo MCA 384 lze používat pouze s předním bezpečnostním panelem.

Otvor dveří je zajištěn pneumatickými pružinami. Operátor musí dveře otevřít zcela, aby byl přístup k přístroji bezpečný/úplný.

Panel ředicí jednotky lze otevřít, aniž by to mělo vliv na provoz zařízení Fluent (s výjimkou zařízení Fluent s instalovaným volitelným světlem UVC – senzor panelu ředicí jednotky při otevření panelu aktivuje rychlé zastavení).

Horní a boční bezpečnostní panely jsou fixní.

2.8.1.1 Přední bezpečnostní panely

Přední bezpečnostní panel zabraňuje přímému přístupu k robotickým ramenům a k prvkům na plošině přístroje během provozu. Slouží účelům osobní bezpečnosti a zlepšuje bezpečnost metod. Přední bezpečnostní panel navíc chrání uživatele před rozlitím vzorku nebo činidla. Existují různé typy předních bezpečnostních panelů.

2 - Bezpečnost Bezpečnostní prvky





Plný přední bezpečnostní panel

Obr. 1: Plný přední bezpečnostní panel

Plný přední bezpečnostní panel má následující funkce:

- Žádný přístup k pohyblivým částem (pohyblivé části, mechanická rizika)
- Ochrana vzorků před vnějšími vlivy (bezpečnost metody)
- Ochrana před rozlitím vzorku nebo činidla



S plnými předními bezpečnostními panely je možné vkládání pouze v dávkách.



Plný přední bezpečnostní panel (UVC)



Obr. 2: Plný přední bezpečnostní panel (UVC)

Plný přední bezpečnostní panel (UVC) má následující funkce:

- Žádný přístup k pohyblivým dílům (pohyblivé díly, mechanická rizika)
- Ochrana vzorků před vnějšími vlivy (bezpečnost metody)
- Ochrana před rozlitím vzorku nebo činidla
- Ochrana před optickým zářením (UVC)



S plnými předními bezpečnostními panely je možné vkládání pouze v dávkách.

A UPOZORNĚNÍ

Pohyblivé díly!

Pohyblivé moduly MCA, FCA a Air FCA mohou způsobit poranění rukou při sahání přes polovinu předního bezpečnostního panelu nebo předního bezpečnostního panelu s expanzí do přístroje za chodu.

Během provozu nesahejte do přístroje.

2 - Bezpečnost Bezpečnostní prvky





Poloviční přední bezpečnostní panel

Obr. 3: Poloviční přední bezpečnostní panel

Poloviční přední bezpečnostní panel má následující funkce:

- Omezený přístup k pohyblivým částem (pohyblivé části, mechanická rizika)
- Ochrana před rozlitím vzorku nebo činidla



S polovičním předním bezpečnostním panelem má obsluha omezený přístup k plošině přístroje. Vkládání a vyjímání podložek je možné bez otevření panelu, tj. operátor má možnost vkládat vzorky nebo činidla během běhu metody.



Přední bezpečnostní panel s rozšířením



Obr. 4: Přední bezpečnostní panel s rozšířením

Přední bezpečnostní panel má následující funkce:

- Omezený přístup k pohyblivým částem (pohyblivé části, mechanická rizika)
- Ochrana před rozlitím vzorku nebo činidla
- Umožňuje použití přední odpadní stanice DiTi, která vyčnívá z plošiny a vyžaduje otvor v předním bezpečnostním panelu směřující dolů.



S předním bezpečnostním panelem s rozšířením je možné vkládání pouze v dávkách.

2.8.1.2 Bezpečnostní panely pro volitelná zařízení

Pokud je na straně produktu Fluent přidáno nebo odstraněno volitelné zařízení, musí být nainstalován příslušný boční bezpečnostní panel. Viz "Zákaznická podpora" [▶ 199].

2.8.2 Zámky dveří přístroje (volitelné)

Dva volitelné zámky dveří slouží k zabránění otevření předního bezpečnostního panelu a chrání probíhající proces. Tím se zabrání nedovolenému přerušení běhu procesu. Chcete-li zastavit proces, můžete zadat požadavek na pozastavení pomocí dotykového displeje.



2.8.3 Zámky dveří skříně

Pokud má dlouhá osa RGA přístup pod plošinu, musí být dveře skříně, které jsou nejblíže místu přístupu, vybaveny volitelným snímačem zámku dveří. Pokud je implementován více než jeden přístupový bod na plošinu, nebo pokud se přístupový bod v průběhu životnosti přístroje změní, musí být každé dveře v blízkosti přístupového bodu vybaveny snímačem zámku dveří.

Pokud je přístroj vybaven digestoří HEPA, musí být všechny dveře skříně vybaveny snímačem zámku dveří.

2.8.4 Optické záření (UVC)

Zařízení Fluent může být vybaveno volitelným odsavačem par HEPA, který obsahuje světlo UVC, nebo samostatným světlem UVC.

Je nutné se vyhýbat záření UVC, protože může vést k poranění. Světlo UVC se automaticky vypne při otevření předního bezpečnostního panelu a v případě volitelného světla UVC také při otevření krytu ředicí jednotky. Ve spojení se světlem UVC se na produkt Fluent instalují speciální bezpečnostní panely odolné proti záření UVC.

Světlo UVC lze použít v postupech dekontaminace. Vhodnost a účinnost použití UVC pro jednotlivé procesy musí ověřit uživatel.



Viz též příručku dodanou výrobcem digestoře HEPA.

2.8.5 Externí zámky dveří

Vnější dveřní zámky budou implementovány u instalací produktu Fluent ve vnějším krytu. Dveřní panely vnějšího krytu nahrazují mechanickou bezpečnostní funkci předního bezpečnostního panelu a dveří skříně produktu Fluent a vnější dveřní doky s integrovanými snímači nahrazují funkci snímače dveří a zámku dveří předního bezpečnostního panelu a dveří skříně produktu Fluent.



Vnější zámky dveří neumožňují funkci ActiveStop. Chcete-li zastavit nebo pozastavit proces, můžete zadat požadavek na pozastavení pomocí dotykového displeje.



2.9 Bezpečnostní značky produktu

K přístroji Fluent jsou z bezpečnostních důvodů upevněny bezpečnostní značky. Poškozené, ztracené nebo nečitelné bezpečnostní značky musí být neprodleně nahrazeny, jak je vidět na obrázku. Význam bezpečnostních symbolů naleznete v části "Konvence bezpečnostních sdělení" [▶ 12].





Obr. 6: Přístroj s prvkem UVC





Obr. 8: Biologické nebezpečí





Přístroj s polovičním předním

panelem

Produkt Fluent je dodáván s bezpečnostní značkou biologického nebezpečí, kterou by měl uživatel použít v případě použití biologicky nebezpečných látek. Štítek nalepte na přední dveře na místo, které je viditelné uživateli a vhodné k použití.



Obr. 9: Přístroj s polovičním předním bezpečnostním panelem

bezpečnostním

2 - Bezpečnost Bezpečnostní značky produktu

Přístroj s předním bezpečnostním

panelem s rozšířením





Obr. 10: Přístroj s předním bezpečnostním panelem s rozšířením

MCA



Obr. 11: Bezpečnostní značka na modulu MCA 384




Obr. 12: Rozšíření plošiny

2.9.1 Pracovní stanice míchání a propichování

Bezpečnostní kryt FCA



Obr. 13: Bezpečnostní kryt



2.10 Laserové záření

Produkt Fluent může být vybaven laserovými skenery čárových kódů. Laserové záření z těchto skenerů čárových kódů je kolimovaný svazek laseru s nízkým výkonem ve viditelném spektru. Třídy laseru každého skeneru čárového kódu a celého systému Fluent jsou uvedeny na bezpečnostním štítku laseru připevněném k příslušnému hardwaru.

Všechny moduly s lasery jsou označeny příslušnými bezpečnostními štítky.

Přístroj Fluent byl testován a certifikován podle IEC 60825-1:2007 a IEC 60825-1:2014.



A UPOZORNĚNÍ

Produkt Fluent je laserový výrobek třídy 1 podle normy IEC 60825-1:2014, který vyzařuje laserové záření.

Laserový paprsek může způsobit oslnění, zábleskovou slepotu a paobrazy.

• Nedívejte se do laserového paprsku ani do jeho odrazů.

2.10.1 Zařízení s laserovým zářením

Na zařízení lze namontovat samostatný skener čárových kódů.

Dbejte na to, aby byl bezpečnostní štítek vždy správně připevněn ke skeneru čárových kódů:

- Vysvětlující štítek laserového záření (A): Označuje LASEROVÝ VÝROBEK TŘÍDY 2 podle normy IEC 60825-1, který obsahuje vestavěný viditelný laserový skener čárových kódů s nízkým výkonem. Dává uživateli pokyn, aby se nedíval do laserového paprsku nebo jeho odrazu.
- Lasery třídy 2 jsou provozovány pouze za chodu systému a nemají žádné rozhraní pro operátora.

Umístění štítku	Vysvětlení
A LASER RADIATION DO NOT STARE INTO BEAM Dutput 8500 Putse duration 1120 CLASS 2 LASER PRODUCT IEC/EN 60825-1 : 2007	Samostatný skener čárových kódů namontovaný na hotelu: štítek umístěný pod skenerem.
LASER RADIATION Do not State into BEAM Output 05,000 Wavelength 65,000 Wavelength 65,000 Utput 05,000 Wavelength 65,000 Utput 05,000 Utput 05,000 Utput 05,000 Wavelength 65,000 Utput 05,000 Utput 05,0	Samostatný skener čárových kódů namontovaný na robotickém upínacím ramenu: štítek umístěný na skeneru.



Umístění štítku	Vysvětlení
	ID vkládání: štítek umístěný na zadní straně skříně skeneru.
	ID vkládání: štítek umístěný na boční straně skříně skeneru.

2.11 Prohlášení o dekontaminaci

Kromě pravidelné péče o systém a v souladu se standardními laboratorními předpisy musí být přístroj Fluent a jeho součásti a příslušenství důkladně dekontaminovány za následujících okolností:

- Před zahájením prací údržby nebo servisu na přístroji Fluent a především před zásahem FSE na přístroji Fluent
- V případě nehod (např. srážka, rozlití apod.)
- Před vrácením přístroje Fluentnebo jeho součástí nebo příslušenství společnosti Tecan (např. za účelem opravy)
- Před uložením
- Před likvidací
- Obecně před přesunem přístroje Fluent nebo jeho částí z umístění

Majitel přístroje nese plnou odpovědnost za účinnou dekontaminaci veškerého zařízení.

Před zásahem FSE na přístroji Fluent a před vrácením přístroje Fluent nebo jeho součástí nebo příslušenství společnosti Tecan musí majitel přístroje vyplnit a podepsat formulář Prohlášení o dekontaminaci a potvrdit, že dekontaminace byla provedena v souladu s pokyny správné laboratorní praxe. Pro získání tohoto formuláře se obraťte na místní servisní organizaci a nahlédněte do části Dekontaminace.



Společnost Tecan si vyhrazuje právo odmítnout zabývat se přístrojem Fluent nebo jeho součástmi či příslušenstvím, ke kterým není přiložen formulář Prohlášení o dekontaminaci.

2.12 Hlášení incidentu

Jakýkoli závažný incident, ke kterému došlo v souvislosti se zařízením, musí být oznámen výrobci a příslušnému orgánu členského státu, kde se uživatel nebo pacient nacházejí. Adresu výrobce naleznete v části Výrobce.



3 Technické údaje

3.1 Typový štítek



- Vor Service oder Wartungsarbeiten Netzstecker ziehen
- Prior to any repair or maintenance job disconnect mains power cord .
- Avant tout type d'intervention, retirer la prise de raccordement au secteur et lire attentivement le manuel
- Prima di esegurie qualsiasi lavoro di manutenzione o servizio, disconnettere il cavo di alimentazione dalla presa di corrente
- Antes de cualquier intervención de servicio o mantenimiento apagar y desconectar el instrumento
- https://www.tecan.com/manuals



(01)07640137481124(11)991231(21)9912123456



VICLAST LASER PRODUCT, THIS PRODUCT COMPLIES WITH 21 CFR 1040.10 AND 1040.11 EXCEPT FOR CONFORMANCE WITH IEC 60825-1 Ed. 3., AS DESCRIBED IN LASER NOTICE No. 56, dated May 8, 2019. IEC 60825-1:2014"



Obr. 14: Typový štítek

Typový štítek se nachází na zadní straně produktu Fluent a obsahuje následující informace:

ldentifikační údaje	Model
	REF: Objednací informace (číslo materiálu a úroveň revize)
	Datum výroby (RRRRMMDD)
	SN: Sériové číslo
Technické údaje	U, f: Napájecí napětí (V), frekvence (Hz)
	P: Příkon (W)
Údaje adresy	Název a adresa výrobce
Údaje o shodě	Označení shody



3.2 Štítek se sériovým číslem



Obr. 15: Štítek se sériovým číslem

Štítek se sériovým číslem je umístěn uvnitř krytu na pravé straně zadní části přístroje a uvádí následující údaje:

ldentifikační údaje	Model
	REF: Objednací informace (číslo materiálu a úroveň revize)
	SN: Sériové číslo
Údaje adresy	Název a adresa výrobce





3.3 Rozměry a hmotnost

	Rozměr	Fluent 480	Fluent 780	Fluent 1080
а	Celková délka	1150 mm (45,28 in)	1650 mm (64,96 in)	2150 mm (84,65 in)
b	Půdorysná hloubka	780 mm (30,71 in)		
с	Celková hloubka	923 mm (36,34 in)		
d	Celková výška na skříni	1977 mm (77,8 in)		

Komponenta	Fluent 480	Fluent 780	Fluent 1080
Základní jednotka	120 kg (264,5 lb)	140 kg (308,6 lb)	190 kg (418,9 lb)
Balení	61 kg (135 lb)	83 kg (183 lb)	106 kg (234 lb)
FCA		10,4 kg (22,9 lb)	
MCA 384	12,6 kg (27,8 lb)		
384kanálová hlava	7,2 kg (15,9 lb)		
RGA	10,2 kg (22,4 lb)		
RGA-Z	10,6 kg (23,4 lb)		



Komponenta	Fluent 480	Fluent 780	Fluent 1080
cXP		1,2 kg (2,6 lb)	
MCA 96 včetně volitelného upínadla		19,7 kg (43,43 lb)	

3.4 Napájení

OZNÁMENÍ

Přehřátí napájecího zdroje

Může dojít k poškození nebo zničení napájecího zdroje.

- Napájecí zdroj nesmí být zakrytý.
- Musí být zajištěn odvod tepla z napájecího zdroje.



K napájení nesmí být připojena externí zařízení. Mohou vést k resetování nebo zastavení systému Fluent

Tab. 2: Plynulý příkon

Napájení	Jmenovitý výkon
Linkové napětí (jedna fáze)	100 až 240 V AC
Vstupní proud	9,8 A (při 100 V) – 4 A (při 240 V)
Frekvence	50–60 Hz

Tab. 3: Plynulý výkon

Napájení	Jmenovitý výkon
Výstupní napětí	24–28 V nastaveno ve výrobě: 25,2 V
Nepřetržitý výkon	500 W
Špičkový výkon (časový limit)	1500 W po dobu 3 sekund
Hmotnost	3,8 kg (8,5 lb)

Maximální kolísání síťového napětí: ±10 % jmenovitého napětí.

Klasifikace z hlediska elektrické bezpečnosti podle norem EN/IEC:

Tab. 4: Elektrické specifikace (bezpečnost)

Kategorie přepětí	II	IEC 60664-1
Stupeň znečištění	2	(EN) IEC 61010-1





3.5 Datová a napájecí připojení

Α С Napájecí jednotka

- 11 Rozhraní USB

D Nástěnná zásuvka

12 Napájecí kabel

Na obrázku jsou znázorněny součásti vzorového systému s datovými a napájecími přípojkami. Součásti přístroje Fluent jsou zobrazeny v obdélníku. Vypínač napájení přístroje je součástí napájecí jednotky. Napájecí kabel se připojuje do síťové zásuvky.

Veškeré datové přenosy do zařízení Fluent a z něj probíhají přes rozhraní USB. Kabel USB je připojen k počítači, který přístroj řídí.

3.6 Podmínky prostředí

▲ UPOZORNĚNÍ

Nesprávné pipetovací objemy

Výsledky pipetování mohou být ovlivněny provozními podmínkami.

Kondenzace může mít vliv na elektronické součástky.

Jestliže je produkt Fluent skladován nebo přepravován při teplotách nižších než je pokojová teplota, bude po instalaci vyžadovat několik hodin aklimatizace.





Provozní podmínky

Podmínky přepravy

Podmínky skladování Produkt Fluent je určen pouze pro provoz a skladování v interiéru.

Provozní teplota	15–32 °C (59–90 °F)
Provozní vlhkost	30–80 % relativní (nekondenzující) při 30 °C (86 °F)
Provozní nadmořská výška	max. 2000 m nad hladinou moře

Provozní podmínky pro manipulaci s kapalinami a pipetování:

Pokojová teplota	20–25 °C (68–77 °F)
Provozní vlhkost	30–60 % relativní (nekondenzující)
Provozní nadmořská výška	kolem 500 m nad hladinou moře
Odpařování	Prostředí se zvýšeným prouděním vzduchu (v důsledku laminárního proudění, klimatizace nebo ventilace atd.) zvyšuje riziko odpařování, které může snížit přesnost pipetování, zejména u malých objemů nebo těkavých látek. OZNÁMENÍ Zajistěte, aby podmínk validace odpovídaly podmínkám provozu.
Přepravní teplota	-20 až 60 °C (-4 až 140 °F)
Přepravní vlhkost	20–80 % relativní (nekondenzující)
Teplota skladování	1–60 °C (34–140 °F)
Skladovací vlhkost	5–80 % relativní (nekondenzující) při

30 °C (86 °F) nebo nižší

3.7 Emise a odolnost

Emise hluku < 60 DBA (akustický tlak), měřeno ve vzdálenosti 1 m od přístroje. Hladina hluku může v průběhu procesu krátkodobě překročit 78 dB.

EMC Produkt Fluent splňuje požadavky na emise a odolnost popsané v IEC 61326-1 a IEC 61326-2-6. Před zahájením provozu produktu Fluent by však mělo být vyhodnoceno elektromagnetické prostředí. Provozovatel je odpovědný za to, aby zajistil kompatibilní elektromagnetické prostředí pro přístroj Fluent, aby Fluent fungoval tak, jak má. Přístroj Fluent je klasifikován jako zařízení SKUPINY 1, TŘÍDY B (CISPR 11).

> Zařízení je určeno k použití v ZÁKLADNÍM ELEKTROMAGNETICKÉM PROSTŘEDÍ (IEC 61326-1) a V PROSTŘEDÍ PROFESIONÁLNÍHO ZDRAVOTNICKÉHO ZAŘÍZENÍ (IEC 61326-2-6). Při použití v PRŮMYSLOVÉM ELEKTROMAGNETICKÉM PROSTŘEDÍ



(IEC 61326-1) a v DOMÁCÍM ZDRAVOTNICKÉM PROSTŘEDÍ (IEC 61326-2-6) je pravděpodobné, že bude fungovat nesprávně.

Pokud existuje podezření, že je výkon ovlivněn elektromagnetickým rušením, lze správnou funkci obnovit zvětšením vzdálenosti mezi zařízením a zdrojem rušení.

Nepoužívejte produkt Fluent v těsné blízkosti zdrojů silného elektromagnetického záření (např. nestíněných úmyslných RF zdrojů), protože mohou narušit správnou funkci.

FCC15 Toto zařízení bylo testováno a vyhovuje limitům pro digitální zařízení třídy B podle části 15 pravidel FCC. Limity jsou navrženy tak, aby poskytovaly přiměřenou ochranu proti škodlivému rušení při instalaci v obytných budovách. Toto zařízení generuje, využívá a může vyzařovat vysokofrekvenční energii, a pokud není instalováno a používáno v souladu s pokyny, může způsobovat škodlivé rušení rádiové komunikace. Neexistuje však žádná záruka, že k rušení nedojde v konkrétní instalaci. Pokud toto zařízení způsobuje škodlivé rušení příjmu rozhlasu nebo televize, což lze zjistit vypnutím a zapnutím zařízení, doporučujeme uživateli, aby se pokusil napravit rušení pomocí jednoho nebo více následujících opatření:

- Změňte orientaci nebo umístění přijímací antény.
- Zvětšete vzdálenost mezi zařízením a přijímačem.
- Připojte zařízení do zásuvky v jiném okruhu, než ke kterému je připojen přijímač.
- Obraťte se na prodejce nebo zkušeného rozhlasového či televizního technika.



4 Popis funkce

Tato kapitola vysvětluje základní funkci Fluent, znázorňuje strukturu a poskytuje funkční popis sestav.

4.1 Přehled

Produkt Fluent se používá pro pipetování pomocí robotických ramen. Robotická ramena mohou nasávat a dávkovat do různých nádob, jako jsou zkumavky se vzorky nebo mikrodesky.

Fluent je k dispozici ve třech různých velikostech:

- Fluent 480
- Fluent 780
- Fluent 1080



Obr. 17: Přehled přístrojů (přístroj se může lišit od vyobrazení)

A Plošina

Robotická ramena

C Doplňky a zařízení



K dispozici je také sada pro ochranu před zemětřesením pro oblasti náchylné k zemětřesení.

В

Další informace naleznete v části "Zákaznická podpora" [> 199].

4.2 Plošina

SegmentyPlošina Fluent, která je vzorkovací zónou přístroje, se skládá ze segmentů.
Segmenty plošiny jsou vyměnitelné součásti plošiny, které mohou mít různé
rozměry a vlastnosti. UPOZORNĚNÍ! Neprovozujte systém bez segmentů
plošiny.

Číslo mřížky Šířka segmentu je vyjádřena v číslech mřížky. Mřížka je široká 25 mm a odpovídá vzdálenosti mezi polohovacími kolíky segmentu.



Čísla mřížky se používají také pro vyjádření umístění segmentů nebo podložek na plošině.

4.2.1 Nosiče



Obr. 18: Plošina Fluent



Nosiče jsou komponenty plošiny určené k uložení laboratorního nebo spotřebního materiálu na plošině.

Podložky jsou nosiče, které se nasouvají a sesouvají na segmentech mřížky a obvykle drží zkumavky se vzorky nebo žlábky s činidly.

Segmenty jsou statické prvky zablokované na plošině. Některé segmenty mají soupravy (segmenty souprav), do kterých se ukládá laboratorní vybavení, jako jsou mikrodesky nebo desky s hlubokými nádobami, nebo spotřební materiál, například boxy DiTi. Některé segmenty mají mřížkové kolíky (mřížkové segmenty) pro vkládání a vykládání podložek.



4.2.2 Podnosy plošiny



Obr. 19: Podnos pracovní plochy

Podnosy pracovní plochy, které jsou umístěny pod dynamickými segmenty pracovní plochy, zachycují rozlitou kapalinu, která se může objevit v oblasti ručního zakládání pracovní plochy. Systém by měl být provozován s co největším počtem podnosů pracovní plochy nainstalovaných pod pracovní plochou, aby byly zachyceny všechny rozlité kapaliny. **UPOZORNĚNÍ! Neprovozujte systém bez podnosů a segmentů pracovní plochy.**

Výřezy v podnosech pracovní plochy pro nářadí a nástroje jsou povoleny pouze v kabinetové verzi.

Vyvýšená místa na každém podnosu pracovní plochy jsou navržena tak, že umožňují přítomnost výřezů bez ovlivnění objemu záchytu. Tyto výřezy umožňují použít odpadní skluzy v plošině nebo jiné integrace zařízení při použití verze montované na kabinet. U přístupových bodů v plošině lze podnosy umístit na polici v kabinetu pod integrovanými zařízeními.



Obr. 20: Podnosy pracovní plochy pod segmenty pracovní plochy

Podnosy pracovní plochy nebudou přítomny tam, kde RGA vyžaduje přístup k zařízení pod pracovní plochou. S přístrojem je dodávána sada podnosů pracovní plochy. Podnosy pracovní plochy lze podle potřeby umýt nebo vyměnit. Viz část "Konec dne" [> 116].



4.2.3 Umístění na segmentu



Obr. 21: Blokovací kolíky a polohovací kolíky

Fluent používá kolíky ke správnému umístění podložek, adaptérů nebo doplňků na segmentu. Podložky jsou navrženy tak, aby se nasazovaly na kolíky. Jejich polohu pak lze zkontrolovat přečtením čísla mřížky na přední straně přístroje. Blokovací kolíky drží podložky ve správné poloze.

A Blokovací kolíky B Polohovací kolíky





4.2.4 Pozice segmentu



Obr. 22: Boční pozice a pozice mřížky

A Boční pozice B

Boční pozice (AB, YZ) lze použít k umístění laboratorního vybavení, se kterým manipuluje RGA.

Pozice mřížky



Při použití FCA nebo MCA není možné pipetovat v bočních pozicích.

Očíslované pozice mřížky (1-n) jsou přístupné pro pipetovací ramena. V konfiguracích s více rameny však nejsou všechny číselné mřížky přístupné všem pipetovacím ramenům. V závislosti na konfiguraci ramena přístroje mohou platit určitá omezení.

4.3 Robotická ramena

Produkt Fluent může být vybaven různými robotickými rameny:

- Flexibilní rameno kanálu (FCA Flexible Channel Arm)
- Vícekanálové rameno (MCA 96 a MCA 384 Multiple Channel Arm)
- Robotické upínací rameno (RGA Robotic Gripper Arm)

Robotická ramena lze vybavit různým příslušenstvím.





4.3.1 Flexibilní rameno kanálu (FCA – Flexible Channel Arm)

Obr. 23: Flexibilní rameno kanálu

6

Pokud není povoleno přenášení, důrazně se doporučuje používat jednorázové špičky s filtry.

Rameno FCA (A) je vybaveno pipetovacími špičkami a může řídit manipulaci s kapalinou až pro 8 samostatných kanálů.

Rameno FCA v konfiguraci s adaptéry DiTi má volitelné upínadlo FCA, který umožňuje určité pohyby laboratorního vybavení – viz "Upínadlo FCA" [> 60].

4.3.1.1 FCA s kapalinovým systémem (FCA s kapalinou)

Rameno FCA s kapalinovým výtlačným systémem je naplněno systémovou kapalinou, která je dodávána injekčními čerpadly. Používá se k pipetování kapalin s různými rozsahy objemu v závislosti na použitých špičkách a velikosti stříkačky. FCA s kapalinou lze nakonfigurovat s pevnými omyvatelnými špičkami nebo s jednorázovými adaptéry špičky.



Společnost Tecan doporučuje používat jako systémovou kapalinu deionizovanou vodu.

4.3.1.2 FCA se vzduchovým systémem (FCA se vzduchem)

Rameno FCA se systémem vytlačování vzduchu se používá k pipetování kapalin pohybem pístu uvnitř pipetovacího kanálu. Rameno Air FCA je nakonfigurováno s jednorázovými adaptéry špiček.





4.3.2 Vícekanálové rameno 384 (MCA 384)

Obr. 24: Vícekanálové rameno



Pokud není povoleno přenášení, důrazně se doporučuje používat jednorázové špičky s filtry.

Rameno MCA (A) je robotické rameno s vícekanálovou pipetovací hlavou. Všechny kanály pipetovací hlavy nasávají a dávkují současně. Pipetovací hlava může vyměňovat adaptéry hlavy. Různé typy adaptérů hlavy umožňují různé formáty pipetování:

- Kompatibilní s portfoliem MCA DiTi
- Pipetování do 125 µl
- Rozsah pipetování až 500 µl s adaptérem pro volitelný/rozšířený objem
- · Bez detekce hladiny kapaliny
- MCA 384 s 384 jednorázovými špičkami
- MCA 384 s 96 jednorázovými špičkami (deska adaptéru)
- MCA 384 s 384 fixními, omyvatelnými špičkami
- MCA 384 s 96 fixními, omyvatelnými špičkami

4.3.3 Vícekanálové rameno 96 (MCA 96)

Modul MCA 96 je robotické rameno s vícekanálovou pipetovací hlavou. Všech 96 kanálů pipetovací hlavy nasává a dávkuje současně. Mezi klíčové vlastnosti ramena MCA 96 patří:





Obr. 25: Hlava/upínadlo MCA 96

- Kompatibilní s vodivým portfoliem FCA-DiTi do 1000 μl (včetně)
- Rozsah pipetování od 1 µl do 1000 µl
- Provádí (kapacitní) detekci hladiny kapaliny (cLLD)
- Provádí částečný odběr DiTi a pipetování s posunem (např. pro ředění) pomocí produktů se špičkami s kombinovaným zásobníkem Tecan)
- Volitelné upínadlo pro jednoduché úkony manipulace s laboratorním vybavením

kompatibilita s funkcí cLLD.

Hlava ramena MCA 96 umožňuje používat jednorázové špičky, aby byla zajištěna

Technologie cLLD je s ramenem MCA 96 ověřena pouze pro použití ve vaničkách.

První špička, která vnikne do kapaliny, aktivuje signál cLLD. Nerovnoměrná výška naplnění (například v ručně naplněných jamkách mikrodesky) by mohla vést k nezamýšlenému zpracování a chybným výsledkům.

6

S ramenem MCA 96 používejte pouze nosiče a laboratorní vybavení kompatibilní s funkcí cLLD. Pro částečný odběr DiTi používejte kompatibilní zásobníky: zásobníky Tecan SLAS. Podrobnosti naleznete v referenční příručce (ref. [4]).

Volitelné upínadlo, které nelze upravit na místě, pro rameno MCA 96 je vhodné pro jednoduchou manipulaci s laboratorním vybavením.



Upínadlo lze použít například pro následující úkony:

- Přesouvání mikrodesek a DWP v režimu na šířku do cílové polohy na pracovním stole
- Práce s naskládanými zásobníky DiTi, včetně vyhazování prázdných vrstev do odpadu
- Manipulace s víčky mikrodesek

Upínadlo ramena MCA je zabudováno jako submodul do hlavy ramena MCA, takže představuje kompaktní a cenově výhodnou alternativu ke specializovanému robotickému upínacímu ramenu (RGA). Upínadlo ramena MCA 96 nedokáže provádět složitější úkony manipulace s laboratorním vybavením, jako jsou:

- Přístup k hotelům
- Manipulace s laboratorním vybavením v orientaci na výšku
- Úkony zahrnující otáčení

Tyto složité úkony nadále vyžadují doplnění systému o specializované robotické upínací rameno (RGA). Upínadlo ramena MCA 96 obsahuje stejnou stanici pro výměnu prstů jako upínadlo ramena RGA.

4.3.4 Robotické upínací rameno (RGA – Robotic Gripper Arm)



Obr. 26: Robotické upínací rameno

Rameno RGA (A) je robotické rameno s upínací hlavou a upínacími prsty. RGA přenáší mikrodesky a další laboratorní vybavení mezi pozicemi na plošině, periferními zařízeními a úložným prostorem pro laboratorní vybavení:

Robotické uchopovací rameno standardní výšky (RGA standard Z) může přistupovat k předmětům umístěným na plošině nebo na spodní plošině.

Vysoké robotické uchopovací rameno (RGA long Z) má přístup k předmětům umístěným na plošině, spodní plošině a níže.

4.3.4.1 Robotická upínací hlava

Rameno RGA lze vybavit dvěma různými variantami robotických upínacích hlav.

Běžná upínací hlava nabízí výběr upínacích prstů, které lze ručně vyměňovat.



Systém výměny prstů (FES) nabízí automatickou výměnu prstů s výběrem sad prstů upínadla. Sady prstů jsou namontovány na dokovací stanici umístěné na standardním segmentu soupravy. Sady prstů jsou automaticky odebírány a umísťovány robotickým ramenem. Typ prstu a výměna prstů jsou sledovány. V rámci jedné metody lze použít libovolný prst nebo všechny prsty.

4.3.5 Příslušenství ramena



Obr. 27: Příslušenství ramena

- A Fixní hroty
- **C** Prsty upínadla

- B Jednorázové špičky
- D Skener čárových kódů pro RGA

4.3.5.1 Fixní hroty



Pokud není povoleno přenášení, důrazně se doporučuje používat jednorázové špičky s filtry.

Omyvatelné, opakovaně použitelné špičky pro aspiraci a dávkování jsou k dispozici pro FCA a MCA.

4.3.5.2 Jednorázové špičky

Špičky jsou k dispozici v podnosech nebo boxech (jednotlivě nebo v soupravách) podle typu. Po aspiraci se špičky vyřadí nebo znovu uloží. Špičky se vyhazují pomocí systému vyhazování jednorázových špiček do odpadního žlabu namontovaného na segmentu plošiny.

4.3.5.3 Prsty upínadla

Pro běžnou hlavu upínadla ramena RGA a upínadlo ramena MCA 96, které jsou vybaveny systémem FES (Finger Exchange System) pro výměnu prstů, jsou dostupné různé typy prstů upínadla.

Excentrické Excentrické upínací prsty přenášejí předměty na mikrodesce v rámci pipetovací oblasti i mimo ni. Předměty na desce jsou uchopovány ze stran. K dispozici jsou dvě varianty:

- Standardní délka prstů pro vkládání mikrodesek do hotelů a zařízení.
- Excentrické dlouhé prsty pro vkládání hlubších zařízení, jako je čtyřdrážka, buněčná deska, sledovaný inkubátor.



Excentrické	
prsty upínadla	
ramena MCA 96	

Centrické

upínací prsty

Excentrické prsty upínadla ramena MCA 96 přemisťují objekty na mikrodeskách v pipetovací oblasti. Tento typ prstů dokáže uchopit objekty na deskách pod hlavou upínadla ramena MCA 96. (Tento typ prstů NENÍ kompatibilní s upínadlem ramena RGA.)

Centrické upínací prsty přenášejí předměty na mikrodesce v rámci pipetovací oblasti a pod ní. Předměty na desce jsou uchopovány shora. Platí pouze pro běžnou hlavu upínadla ramena RGA.

Zkumavkové
prstyZkumavkové prsty přenášejí předměty ve zkumavkách v rámci pipetovací oblasti
a pod ní. Platí pouze pro běžnou hlavu upínadla ramena RGA.

4.3.5.4 Skener čárových kódů

Rameno RGA může být vybaveno horizontálním skenerem čárových kódů na mikrodeskách a boxech DiTi.



Bezpečnostní pokyny pro třídu laseru je potřeba pečlivě si přečíst a dodržovat je. Viz též příručku dodanou výrobcem skeneru čárových kódů.

4.4 Kapalinový systém (FCA s kapalinovým systémem)



Obr. 28: Kapalinový systém (FCA s kapalinovým systémem)

Kapalinový systém je navržen pro účinné vnitřní i vnější mytí pevných pipetovacích špiček.



4.5 Mycí systém (MCA 384)



Obr. 29: Mycí systém (MCA)

Mycí blok (A) nainstalovaný na segmentu ramena MCA 384 promývá špičky adaptéru fixních špiček po každém pipetovacím cyklu.

4.6 Doplňky a zařízení



Seznamu příkladů doplňků a zařízení třetích stran pro integraci do systému Fluent. Některé doplňky společnosti Tecan a zařízení třetích stran, která lze používat s přístrojem Fluent, jsou určena jen pro výzkumné použití (RUO).

V této části jsou doplňky a zařízení určené pouze pro výzkum označeny hvězdičkou (*).

Další informace naleznete v části "Zamýšlené použití" [▶ 8].

Pasivní doplňky

Aktivní doplňky

- Hotel (úložné zařízení desek)
- Skříň
- Prachový kryt
- Upínadlo FCA
- HEPA digestoř
- Stohovač Fluent
- MIO2
- Te-Shake
- Te-VacS
- Karusel Fluent
- Věž mycího a doplňovacího centra (WRC Wash and refill center) (např. mycí stanice MCA)
- Resolvex i300
- Propichovací špičky a rotátor zkumavek
- Čtečka FRIDA

Další informace naleznete v části "Referenční dokumenty" [▶ 9].

Myčky založené na systému HydroControl



- Zůstatky založené na standardu MT-SICS úrovně 1
- Zařízení kompatibilní se SiLA*
- Agilent Sealer*
- Inheco ODTC

.

- Inheco Heating Cooling s využitím řídicí jednotky MTC/STC
- Cytomat 10*, 20*, 200* a 6000*

Čtečky čárových kódů

Čtečky čárových kódů řady Keynence BL-1300

Skener čárových kódů zkumavek Fluent ID

Čtečky

- Čtečky Tecan řízené systémem Magellan
- Spark a SparkControl Magellan*
- Stolní čtečka Ziath 2D*



Nahlédněte také do příruček dodaných výrobcem doplňku, přístroje nebo zařízení třetí strany. Pokyny je potřeba pečlivě si přečíst a dodržovat.

4.6.1 Skener čárových kódů zkumavek Fluent ID



Obr. 30: Fluent ID

- A Kryt skeneru
- C Nakládací prostor
- E Reflektor

- B Las
 - Laserový skener čárových kódů
 - LED diody

Fluent ID je volitelný modul, který lze použít ke snímání štítků s čárovým kódem zkumavek při vkládání zkumavek na plošinu. Každý modul Fluent ID obsahuje šest vyhrazených mřížkových pozic pro nakládání a snímání štítků s čárovým kódem až šesti podložek. Reflektor se používá k detekci prázdných pozic trubek v podložce. Grafické rozhraní na dotykovém monitoru poskytuje pokyny pro ovládání systému Fluent ID.

D

Laserové záření z tohoto skeneru čárových kódů je kolimovaný svazek laseru s nízkým výkonem ve viditelném spektru s následujícími vlastnostmi.



- Vlnová délka: 655 nm
- Trvání pulzu: 150 µs
- Maximální výkon výstupní energie: 1,0 mW

4.6.1.1 Podložky pro kompatibilní zkumavky Fluent ID

Podložky zkumavek Fluent ID jsou určeny vždy pro jeden typ zkumavek:

- Podložka s 32 pozicemi pro zkumavky s průměrem 10 mm
- Podložka s 32 pozicemi pro zkumavky s průměrem 13 mm
- Podložka s 26 pozicemi pro zkumavky s průměrem 16 mm
- Podložka s 32 pozicemi pro zkumavky 2ml Eppendorf Safe-Lock



Volitelnými zátkami lze zablokovat dvě pozice 26polohové podložky, aby bylo možné ji použít jako 24polohovou podložku, což umožňuje paralelní pipetování ze zkumavek v násobcích osmi.

4.6.2 Upínadlo FCA

Přehled



Obr. 31: Upínadlo FCA

Upínadlo FCA je volitelný doplněk pro rameno FCA nakonfigurované s adaptéry DiTi, které kromě pipetování umožňuje FCA provádět některé pohyby s laboratorním vybavením. FCA může automaticky získat a upustit prsty upínadla FCA za chodu.



Prsty upínadla FCA



Obr. 32: Prsty upínadla FCA

Prsty upínadla FCA je potřeba vyměnit po 2 letech nebo 20000 cyklech použití (jeden cyklus je definován jako vyzvednutí, použití a zaparkování). Cykly se sledují pomocí čítače definovaného v softwaru Fluent Control.

Souprava dokovací stanice upínadla FCA



Obr. 33: Souprava dokovací stanice upínadla FCA

Souprava dokovací stanice upínadla FCA slouží k uložení prstů upínadla FCA. Lze ji namontovat jako standardní soupravu pro mikrodesky na segment plošiny.



4.6.3 Míchání a propichování

Pracovní stanice pro míchání a propichování Fluent je určena pro aplikace, při nichž se kapalina přenáší ze zkumavek na vzorky uzavřené gumovým uzávěrem do zkumavek s gumovým uzávěrem, kdy není nutné uzávěr odstranit a stačí jej propíchnout.



Uzávěr zkumavky je možné propíchnout pouze jednou. Opakované propíchnutí stejné zkumavky není možné.

Propichování bylo testováno u zkumavek BD Vacutainer® s uzávěrem Hemogard a u zkumavek Greiner Vacuette® s hladkým stahovacím uzávěrem a bezpečnostním šroubovacím uzávěrem. Propichování u zkumavek s gumovými zátkami/uzávěry není možné.



Propichovací špičky je nutno pravidelně vyměňovat. Použití bylo ověřeno při počtu 20 000 propíchnutí na jednu špičku.

Pracovní stanice pro míchání a propichování Fluent je v závislosti na velikosti základní jednotky Fluent konfigurována až se 2 kapalinovými rameny FCA, stanicí pro hloubkové mytí a až 4 rotátory zkumavek. Rotátory zkumavek lze integrovat na jakoukoli velikost základní jednotky Fluent a podporují čtení čárových kódů zkumavek, míchání vzorků, propichování a alikvotaci. Další informace o podporovaných typech zkumavek naleznete v části "Podložky rotátoru zkumavek" [» 63].

Pracovní postup lze rozdělit do následujících kroků:

- 1. Čtení čárových kódů při vkládání zkumavek
- 2. Míchání obsahu zkumavky
- 3. Propichování a manipulace s kapalinou pomocí ramena FCA v rotátoru zkumavek s propichovacími špičkami
- 4. Mytí a dekontaminace propichovacích špiček v hloubkové mycí stanici a dekontaminačních žlabech
- 5. Opakování kroku 2 a pokračování
- 6. Na konci skriptu se doporučuje provést prodloužené proplachování mycí stanice včetně centrálního odpadu a předního a zadního čističe.

4.6.3.1 Rotátor zkumavek

Hlavním účelem modulu rotátoru zkumavek je promíchat kapalný obsah zkumavek a posloužit jako nosič pro úkony propichování a pipetování. Jeden rotátor zkumavek má kapacitu 5 podložek rotátorů zkumavek, každý s 24 zkumavkami (tj. celková kapacita 120 zkumavek).

Zařízení obsahuje následující dílčí součásti:

- Integrovaný skener čárových kódů na zkumavkách pro snímání čárových kódů vzorků během vkládání
- Volitelná hluboká mycí stanice s hlubokými žlábky pro dekontaminaci propichovacích hrotů a držák chybových zkumavek. Držák chybových zkumavek lze použít k uložení vzorků v případě chyb při propichování. Mycí stanice je umístěna vedle rotujícího bubnu.



- Rotační buben s dolním držákem zkumavek, který pojme až pět podložek rotátoru zkumavek. Buben provádí míchání vzorku buď rotací o 360°, nebo oscilací s různými úhly a rychlostmi. Držák (kryt) podporuje proces propíchnutí.
- Rotátor zkumavek s rotujícími zkumavkami podporuje kapacitní detekci hladiny kapaliny před a po aspiraci a také po dávkování kapaliny (kontrola přítoku kapaliny) přes uzavřené zkumavky (volitelné nastavení).
- Rotátor zkumavek nainstaluje FSE a hlavní operátor ani uživatel s ním nesmí hýbat.

4.6.3.2 Podložky rotátoru zkumavek

Podložky rotátoru zkumavek jsou určeny k použití na rotátoru zkumavek a k podpoře funkce propíchnutí. Pro podporované typy zkumavek k propíchnutí jsou k dispozici různé podložky zkumavek:

- Podložka rotátoru zkumavek BD 13 × 75 mm, 24 pozic pro zkumavky
- Podložka rotátoru zkumavek BD 13 × 100 mm, 24 pozic pro zkumavky
- Podložka rotátoru zkumavek Greiner 13 × 75 mm, 24 pozic pro zkumavky
- Podložka rotátoru zkumavek Greiner 13 × 100 mm, 24 pozic pro zkumavky
- Podložka rotátoru zkumavek 16 × 100 mm, 24 pozic pro zkumavky



Vícenásobná aplikace je podporována pouze u zkumavek Greiner Vacuette® s hladkým stahovacím uzávěrem a bezpečnostním šroubovacím uzávěrem.

Tab. 5: Kompatibilita zkumavky a podložky

Výrobní linka	Zkumavka			Podložka	
	Průměr [mm]	Délka [mm]	Definice laboratorního vybavení	Kompatibilní podložka	Barva mostu
Greiner Vacuette	13	100	13 × 100 mm Greiner Vacuette s přepážkou	1 × 24 13 × 100 mm podložka rotátoru zkumavek Greiner	šedá
	13	75	13 × 75 mm Greiner Vacuette s přepážkou	1 × 24 13 × 75 mm podložka rotátoru zkumavek Greiner	
	16	100	16 × 100 mm Greiner Vacuette s přepážkou	1 × 24 16 × 100 mm podložka rotátoru zkumavek	černá
BD Vacutainer	13	100	13 × 100 mm BD Vacutainer s přepážkou	1 × 24 13 × 100mm podložka rotátoru zkumavek BD	bílá
	13	75	13 × 75 mm BD Vacutainer s přepážkou	1 × 24 13 × 75 mm podložka rotátoru zkumavek BD	
	16	100	16 × 100 mm BD Vacutainer s přepážkou	1 × 24 16 × 100 mm podložka rotátoru zkumavek	černá



4.6.3.3 Ochrana propichovacího hrotu

Obr. 34: Ochrana propichovacího hrotu

Ochrana propichovacího hrotu je krytka, která slouží k zakrytí ostrého hrotu pro propichování při výměně hrotu a řešení problémů. Chrání uživatele před zraněním a hroty před poškozením.



Ochrana propichovacího hrotu je určena pouze pro jednorázové použití. Po použití musí být všechny ochrany propichovacího hroty vyhozeny do kontejneru na biologický odpad.





4.6.3.4 Nástroj pro odstranění propichovacího hrotu

Obr. 35: Nástroj pro odstranění propichovacího hrotu

Nástroj pro odstranění propichovacího hrotu slouží k zatáhnutí propichovacího hrotu zaseknutého ve zkumavce, který nelze zatáhnout pomocí softwarových příkazů.

4.6.4 Čtečka Frida



Obr. 36: Čtečka Frida



Čtečka Frida je určena k automatické kvantifikaci a normalizaci nukleových kyselin. Vzorky, které se mají měřit pomocí čtečky Frida, je třeba zchladit na 4 °C, aby odpařování vzorku nezhoršilo výsledky měření.

▲ UPOZORNĚNÍ

Vibrace mohou způsobit chybné výsledky!

Vibrace kapky vzorku mohou být příčinou nesprávných výsledků měření a zhoršení bezpečnosti nebo klinického stavu vzorku pacienta.

- Předpokladem vhodného místa instalace je stabilní podlaha.
- Během měření pomocí čtečky Frida nesmí být v blízkosti žádné vnitřní ani vnější zdroje vibrací.
- Vyhněte se zdrojům s rezonanční frekvencí. Zejména je třeba se vyhnout vibracím kolem 36 Hz (2160 ot./min.) a kolem 42 Hz (2520 ot./min.), protože se jedná o rezonanční frekvence zavěšené kapky.

A UPOZORNĚNÍ

Osvětlení místnosti může být příčinou nesprávných výsledků!

Osvětlení místnosti nad modulem může rušit měření, být příčinou nesprávných výsledků měření a zhoršení bezpečnosti nebo klinického stavu vzorku pacienta.

 Robotický systém musí mít neprůhledný horní kryt, přední a zadní panel, aby bylo odstíněno okolní světlo v místě měření pomocí čtečky Frida.

4.6.5 Separátor fází

Separátor fází je určen k detekci separačních fází mezi kapalinami různé viskozity. Jako takový je nezávislý na tom, zda je separační fáze viditelná vně laboratorního vybavení. Separátor fází lze použít v aplikacích, které vyžadují čistý přenos kapalné fáze ze zdrojového do cílového laboratorního vybavení.

Následující příklad představuje typický pracovní postup:

- Odstředění zdrojového laboratorního vybavení s kapalnou směsí za účelem vytvoření odlišné fáze mezi kapalinami.
- Kapaliny se musí lišit viskozitou, aby se během odstřeďování mohla vytvořit fáze.
- Umístěte odvíčkované zkumavky na plošinu zařízení Fluent a k zajištění plné sledovatelnosti použijte čtečku čárových kódů (např. Fluent ID). Dbejte na to, abyste při umisťování nenarušili vrstvu mezi fázemi / kapalnými frakcemi.
- Spusťte protokol definovaný pro separaci frakcí. Funkce separátoru fází modulu Air FCA rozpozná fázi mezi kapalinami a zahájí přesun sledované frakce do cílového laboratorního vybavení.

Ze zdrojového laboratorního vybavení lze extrahovat více než jednu fázi. Je nutné odebrat alespoň část horní fáze, aby nedošlo k rozlití kapaliny (přetečení laboratorního vybavení) na plošinu a k potenciální kontaminaci pipetovacího kanálu nad jednorázovou špičkou během detekce fází.

Technické údaje jsou uvedeny v referenční příručce. Podrobnosti o softwaru FluentControl naleznete v příručce k aplikačnímu softwaru. Viz "Referenční dokumenty" [▶ 9].



Kapalné fáze musí být zřetelně odděleny. Pro separaci plné krve na plazmu a krvinky je důležitá kvalita a předběžná úprava vzorků. Mezi parametry, které mohou ovlivnit detekci fází ve vzorcích krve, patří kvalita vzorku (lipémie, hemolýza), doba skladování, teplota při skladování, podmínky při přepravě, podmínky při odstřeďování (čas, rcf, teplota, náběh, typ rotoru), deformace fáze po odstředění atd.

Pro dosažení optimálních výsledků separace fází by měly být krevní vzorky zpracovány co nejrychleji po odběru. Při zpracování a skladování vzorků je nutné dodržovat podmínky doporučené výrobcem příslušných zkumavek.

Odstřeďování při 2500 rcf po dobu 10 minut při pokojové teplotě se zpomaleným náběhem vede k čisté separaci fází u vzorků plazmy (podmínky interního testování společnosti Tecan).

Zkumavky by neměly být naplněny až po okraj, aby nedocházelo k rozlití kapaliny ze zkumavek během aspirace a detekce fází. Detekce fází obvykle vyžaduje rychlý pohyb ve zkumavce směrem dolů v kombinaci s nízkou rychlostí aspirace, což vede ke zvýšení hladiny kapaliny během detekce.

U zkumavek s vysokou úrovní naplnění je doporučeno před zahájením detekce fází odebrat kapalinu z horní strany.

4.6.6 Resolvex i300



Obr. 37: Modul pracovní desky Resolvex i300



Resolvex i300 je modul, který je možné integrovat do platformy přístroje Fluent. Platforma přístroje Fluent pak může manipulovat s kapalinami a přenášet laboratorní vybavení z pracovní desky přímo na modul Resolvex i300, a to vše v rámci jediného softwarového prostředí.

Resolvex i300 je laboratorní přístroj, který umožňuje automatizovat celou řadu laboratorních procesů pomocí následujících hlavních funkcí:

- Tlakování filtračních kolon (plně nebo částečně naplněných) vzduchem nebo dusíkem
- Stohování a rozkládání filtračních desek na sběrných deskách s integrovaným zdvihacím zařízením pro laboratorní vybavení
- Ochrana vzorků před křížovou kontaminací při stohování a rozkládání filtračních a jímacích prvků pomocí integrované ochrany proti odkapávání
- Dávkování kapalin do filtrů nebo do jímacích prvků (volitelně)
- Odpařování kapalin z jímacích prvků s ohřátým plynem, například vzduchem nebo dusíkem (volitelně)

Mezi typické pracovní postupy při použití přístroje Resolvex i300 patří následující laboratorní procesy:

- Extrakce pevné fáze (SPE) v pracovních postupech přípravy vzorků s použitím hmotnostní spektrometrie
- Podporovaná extrakce kapalin (SLE)
- Další pracovní postupy filtrace, čištění nebo koncentrace využívající přetlak

Další údaje a informace o veškerých postupech s přístrojem Resolvex i300 naleznete v návodu k obsluze k přístroji Resolvex i300. Viz "Referenční dokumenty" [▶ 9].



5 Řídicí prvky

5.1 Ovládací prvky



Dotyková obrazovka zobrazuje metody a popisy a umožňuje operátorovi ovládat přístroj.



5.2 Uživatelské rozhraní



- A Navigační cesta B Pracovní oblast
- C Tlačítka zobrazení / možností / akcí

V uživatelském rozhraní FluentControl má operátor přístup ke spouštění metod obsluhy a péče o systém.

5.2.1 Navigační cesta

Navigační cesta slouží k pochopení hierarchie FluentControl a k navigaci v této hierarchii.

Tlačítko	Název	Funkce
₫ >	Domů	Stisknutím tlačítka se vrátíte na domovskou stránku.
System Care 🛞	Navigační okno	Zobrazení aktuálních a předchozích voleb.
<	Rozšíření nabídky	Stisknutím tlačítka pro rozšíření nabídky zobrazíte možnosti, jako je ovládání světel a přepínání operátorů.

Tab. 6: Tlačítka navigační cesty



5.2.2 Pracovní oblast

Přístup k metodám a popisům pomocí pracovní oblasti uživatelského rozhraní. Zde se zobrazují také podrobnosti o stavu běhu metody.

Tab. 7: Tlačítka pracovní oblasti

Tlačítko	Název	Funkce
	Běh	Stisknutím tlačítka spustíte vybranou metodu.
+	Přidat	Stisknutím tlačítka přidáte do seznamu rychlého spuštění další metody.
Assay 1 unknown	Vybraná metoda	Aktuálně vybraná metoda, která se spustí po stisknutí tlačítka Pokračovat.
Assay 3 Unknown	Dostupná metoda	Metoda, kterou lze vybrat kliknutím.
Start of Day Due: is not executed	Tlačítko rychlého spuštění	Stisknutím tlačítka okamžitě spustíte vybranou metodu.

Tab. 8: Displej pracovní oblasti

Displej	Funkce displeje
00:16:52 10:09 10:43 Vot Time Tot Time	Zobrazuje stav a zbývající čas pro běh metody.
Assay 1 is ready to be started.	Popis aktuálně vybrané metody nebo další informace o aktuální akci.

5.2.3 Tlačítka zobrazení, možností a akcí

Tab. 9: Tlačítka zobrazení, možností a akcí

Tlačítko	Název	Funkce
✓ Ok	Ok	Stiskněte pro potvrzení.



Tlačítko	Název	Funkce
× Cancel	Storno	Stiskněte pro stornování.
Continue	Pokračovat	Stiskněte pro pokračování.
II Pause	Pauza	Stisknutím vyžádejte pozastavení běhu na konci aktuální akce.
Stop	Stop	Stisknutím okamžitě zastavíte běh, a to i uprostřed probíhající akce. Pokud je to možné, systém nabídne možnost obnovení nebo pokračování běhu.
THE Remove	Odebrat	Stisknutím tlačítka odeberete metodu ze zobrazení rychlého spuštění.
View Mode	Režim zobrazení	Stisknutím přepínáte mezi zobrazením seznamu a rychlého spouštění.
Sort by	Seřadit podle	Stisknutím přepnete zobrazení běhu metody mezi abecedním a nejnovějším řazením.

5.2.4 Tlačítka obnovení metody

Tab. 10: Tlačítka zobrazení, možností a akcí

Tlačítko	Název	Funkce
Discard	Vyřadit	Stisknutím vyřadíte stav obnovené metody.
\rightarrow	_	Stisknutím přejdete na další obrazovku.


Tlačítko	Název	Funkce
Recovery Point	Bod obnovení	Stisknutím se vrátíte na předchozí obrazovku ("Bod obnovení").
Run Recovery	Obnovení běhu	Stisknutím pokračujte v běhu.

5.2.5 Tlačítka DeckCheck

Tab. 11: Tlačítka DeckCheck

Tlačítko	Název	Funkce
Left	Levá kamera	Zobrazí snímek pořízený levou kamerou (pouze Fluent 780/1080). Pokud byla u této kamery zjištěna nesrovnalost v uspořádání, zobrazí se na ikoně vykřičník.
Center	Prostřední kamera	Zobrazí snímek pořízený středovou přehledovou kamerou. Pokud byla u této kamery zjištěna nesrovnalost v uspořádání, zobrazí se na ikoně vykřičník.
Right	Pravá kamera	Zobrazí snímek pořízený pravou kamerou (pouze Fluent 780/1080). Pokud byla u této kamery zjištěna nesrovnalost v uspořádání, zobrazí se na ikoně vykřičník.
Pause Alternate	Pozastavení střídání	Na obrazovce se střídají referenční a živé obrázky: Stisknutím tohoto tlačítka při zobrazení referenčního nebo živého obrazu podržíte tento obraz statický.
Resume Alternate	Obnovení střídání	Obraz je statický: Stisknutím tlačítka obnovíte střídání referenčních a živých snímků.



Tlačítko	Název	Funkce
Check	Kontrola	Aktivuje opakovanou kontrolu systému – například když byly provedeny opravy. Zobrazí se výzva k zavření dveří. U tříramenného systému se musí pohybovat prostřední rameno: Pokud nejsou dveře zavřené, kontrola se provede, ale prostřední rameno zablokuje jednu kameru.
Ignore & Continue	Ignorovat a pokračovat	Objeví se pouze v případě, že je tento příkaz v metodě nakonfigurován. Umožňuje ignorovat zvýrazněné nesrovnalosti a pokračovat v provádění skriptu.
Continue	Pokračovat	Objeví se, pokud byly vyřešeny všechny nesrovnalosti nebo pokud systém nenašel žádné nesrovnalosti a byla vybrána možnost vždy zobrazit pro příkaz. Díky tomu mohou být okem viditelné jemné barevné změny, které systém nerozpoznal.

Tab. 12: Displeje

Displej	Popis	Funkce
	Referenční obraz	Referenční obraz je uložen v příkazu skriptu, který zobrazuje požadované rozložení plošiny.
LIVE	Živý obraz	Živý obraz pořizovaný kamerami při běhu scénáře.
	Nesrovnalost (rozdíl oproti referenčnímu obrázku)	Červenými čtverci jsou označeny oblasti, kde byly zjištěny rozdíly mezi referenčními a živými obrazy. Označená oblast může obsahovat více než jednu chybu.





5.3 Chybové signály a stav přístroje



Stavové kontrolky indikují stav přístroje pomocí různých barev, stálých nebo blikajících světel. Horní stavová kontrolka je aktivní pouze při spuštěném softwaru.

Tab. 13: Světelné signály stavových kontrolek

Signál	Barva	Režim	Stav přístroje
	_	vур	Přístroj je vypnutý (odpojený od napájení).
	bílá	"srdeční tep"	Přístroj je zapnutý (řídicí software je připojen, moduly ještě nejsou inicializovány).
	bílá (jen kontrolka napájení)	kontinuální	Stav "zapnutí" přístroje (řídicí software není připojen).



Signál	Barva	Režim	Stav přístroje
	barevné schéma uživatelského rozhraní FluentControl	"srdeční tep"	 Režim nečinnosti Všechny moduly jsou inicializovány; přístroj je připraven ke spuštění metody. Přibližně po jedné hodině v režimu nečinnosti se přístroj přepne do pohotovostního režimu. Pohotovostní režim Všechny osy jsou zabrzděné. Ramena nejsou v ZeroG a nelze s nimi ručně pohybovat. Chcete-li přístroj aktivovat, spusťte metodu nebo požádejte hlavního operátora o výběr nástroje pohybu pro režim ZeroG.
	žlutá	kontinuální	Režim učení Přístroj se "učí" pozice. V tomto režimu může uživatel pohybovat robotickými rameny ručně.
	zelená	kontinuální	Je spuštěna metoda (skript nebo proces). Toto je běžný " výrobní" režim .
	červená	blikající	Chybový stav Na obrazovce řídicího počítače nebo na dotykovém displeji se zobrazuje chybové hlášení.
	barva je uživatelsky konfigurovatel ná	blikající	Výzva pro uživatele Systém čeká na interakci uživatele.
	zelená	blikající	Aktivní zastavení Toto je úmyslná pauza vyvolaná řídicí jednotkou runtime nebo otevřením bezpečnostního panelu. Přístroj se pozastaví, aby umožnil interakci uživatele s plošinou. Operátor může obnovit metodu.





5.4 Stavové LED kontrolky Fluent ID

Obr. 41: LED kontrolky Fluent ID

LED kontrolky Fluent ID signalizují následující stavy:

Tab. 14: LED kontrolky Fluent ID

Signál	Barva	Režim	Stav přístroje
	_	vур	Fluent ID je v nečinnosti.
	bílá	kontinuální	Zapnuté napájení Fluent ID (ale zatím bez inicializace).
	modrá nebo vlastní barva	blikající	Připraveno ke vkládání nebo vykládání podložek.
	zelená	kontinuální	Čárové kódy byly úspěšně naskenovány. Podložka pod dohledem. Nevykládejte, protože tím přerušíte běh.
	červená	blikající	Chybový stav Na dotykové obrazovce se zobrazuje chybové hlášení a požadovaná akce.



6 Operace

6.1 Bezpečnostní pokyny pro tuto kapitolu

A UPOZORNĚNÍ

Chybné výsledky nebo kontaminace přístroje!

Pokud nebyla provedena kvalifikace instalace a kvalifikace provozu nebo pokud nebyly dodrženy provozní postupy uvedené v této příručce, může dojít k nesprávným výsledkům nebo ke kontaminaci přístroje.

- Záznamy o kvalifikaci instalace a provozní kvalifikaci jsou dostupné a známé.
- Metody a postupy, včetně parametrů pipetování, musí být validovány hlavním operátorem.
- Detekce hladiny kapaliny ve spojení s aplikacemi propíchnutí pro FCA a Air FCA musí být ověřena hlavním operátorem.
- Při práci s plnou krví s použitím propichovacích špiček a rotátoru zkumavek a koncentrovaných korozivních dekontaminačních roztoků se doporučuje vyměnit každé tři měsíce obě přípojky mycí stanice (přímo na mycí stanici a v nádobě na systémovou kapalinu), aby nedošlo k jejich ucpání nebo poškození. Potenciálně korozivní tekutiny je nutno před vypouštěním neutralizovat; případně je možné mycí stanici propláchnout neutrální kapalinou (např. vodou).
- Operátor musí být proškolen o provozních postupech, metodách a procesech.



DEPOZORNĚNÍ

Biologická a chemická kontaminace uživatele!

Poškozené prsty upínadla FCA mohou upustit desky. Upuštěné desky mohou způsobit kontaminaci nebezpečnými látkami.

• Po nárazu zkontrolujte prsty upínadla FCA.

A UPOZORNĚNÍ

Ostré hrany a hroty!

Propichovací špičky pracovní stanice míchání a propichování Fluent mají špičaté hroty a ostré hrany, které mohou způsobit poranění.

- Při nakládání přístroje přesuňte rameno FCA do pozice pro uložení pomocí softwarového příkazu.
- Po chybě zakryjte propichovací špičky ochranou propichovacích špiček a ručně přesuňte rameno FCA do polohy uložení. Viz část "Ochrana propichovacího hrotu" [▶ 64].

A UPOZORNĚNÍ

Biologická kontaminace systému!

Na pracovní stanici míchání a propichování Fluent může krev kontaminovat víčka zkumavek.

- Zacházejte se zkumavkami opatrně.
- Používejte ochranné pomůcky.

OZNÁMENÍ

Poškození způsobené korozivními kapalinami!

Vypouštění koncentrovaných korozivních kapalin, jako je např. 2% bělidlo, skrze mycí stanici a hadičky, může způsobit poškození přípojek mycí stanice.

Potenciálně korozivní tekutiny je nutno před vypouštěním neutralizovat; případně je možné mycí stanici propláchnout neutrální kapalinou (např. vodou).

6.2 Provozní režimy

Přístroj Fluent lze spustit ve třech různých provozních režimech:

 Operátor
 Rutinní provozní režim

 • Normální provozní režim, ve kterém se spouští aplikace nebo běžné úlohy péče o systém.

 • Přístroj Fluent je monitorován řídicí jednotkou runtime softwaru FluentControl.

 Hlavní operátor
 Režim definice metody

Tento provozní režim se používá k provádění speciálních úkolů, jako je seřízení za účelem nastavení metody.



FSE

Servisní režim

- Tento provozní režim se používá k provádění speciálních úkolů, jako jsou testy pro zajištění provozní připravenosti.
- Ke spuštění tohoto režimu je vyžadován servisní certifikát.



▲ UPOZORNĚNÍ

Rameno naráží do objektů na pracovním stole

S rameny Fluent lze pohybovat manuálně. Manuální pohyb ramen musí být prováděn jemně, rameno je nutné držet. Nenarážejte rameny do pevných předmětů, a to ani do mechanické zarážky rozsahu ramen

6.3 Uvedení do provozu

6.3.1 Zapnutí přístroje

Chcete-li přístroj zapnout, postupujte následovně:

1. Zapněte napájení vypínačem (A) na zadní straně externího zdroje napájení.



Na zapnutém přístroji bude kontrolka napájení svítit modře. Viz část . Pokud se stavová kontrolka nerozsvítí, spusťte počítač nebo se obraťte na hlavního operátora.

2. Spusťte software FluentControl. Viz část "Spuštění FluentControl" [> 81].



6.3.2 Spuštění FluentControl

- ✓ Provozní postupy musí být dostupné a známé.
- ✓ Záznamy o kvalifikaci instalace a provozní kvalifikaci jsou dostupné a známé.
- ✓ Byla provedena péče o systém.
- ✓ Přístroj je zapnutý.
- Spustte software pomocí Start > Všechny programy > Tecan > FluentControl.

Po několika sekundách se objeví obrazovka Start .

6.3.3 Přihlášení uživatele



Správa uživatelů je k dispozici pouze společně se softwarem Fluent Gx Assurance.

Chcete-li se přihlásit do systému FluentControl, postupujte takto:

- ✓ Software Fluent Gx Assurance je nainstalovaný.
- ✓ Aplikace FluentControl je spuštěná.
- ✓ Správa uživatelů je v aplikaci FluentControl aktivovaná a proces je definovaný.
- 1. Vyberte přiřazený uživatelský profil (A).



2. Zadejte heslo na klávesnici (B).



3. Stiskněte OK (C).



Po přihlášení se přístroj automaticky inicializuje.

6.3.4 Umístění segmentů

Postup umístění segmentů:

- Všechny segmenty, nosiče, doplňky a zařízení musí být umístěny v souladu se zvolenou metodou.
- ✓ Segmenty jsou vyčištěné a v perfektním stavu.
- ✓ Segmenty jsou umístěny na odpovídající pozici mřížky.
- 1. Při instalaci aktivního nosiče modulu MCA 384 připojte kabel.



2. Spusťte segment na zadní část plošiny.



- 3. Vyrovnejte zadní hranu s krytem zadního kanálu nebo s nástavcem přístroje.
- 4. Opatrně spusťte dolů přední část segmentu (A).



 Otočte upínací páku zleva doprava do zavřené polohy. Viz část "Kontrola segmentu" [▶ 192].

6.3.5 Odebírání segmentů

6

Segmenty Fluent ID nejsou navrženy k odebrání! Jsou připojeny přímo k elektronice přístroje. Připojení k elektronice může provádět pouze kvalifikovaný FSE.

Viz část Zákaznická podpora.

Chcete-li odebrat segmenty, postupujte takto:

- Všechna činidla, vzorky, stojany, podložky a desky byly ze segmentu odebrány.
- Na segmentu není nic umístěno.
- 1. Otočte upínací páku (A) zprava doleva do otevřené polohy.







Segment je odemknutý a žlutá značka na upínací páce je vidět.

2. Posuňte segment přibližně o 4 mm dopředu.



3. Zvedněte segment (B) vpředu.



Před odebráním aktivního nosiče MCA odpojte kabel.



4. Segment uložte na čistém a suchém místě, aby nedošlo k jeho poškození.

6.3.6 Vkládání standardních podložek

OZNÁMENÍ

Poškození v důsledku nesprávného vkládání nebo vykládání

Poškození podložek a kolíků.

- Vyrovnejte podložku horizontálně s plošinou.
- Jednou rukou podepřete přední část podložky.
- Při vykládání se ujistěte, že se podložka před zvednutím uvolnila ze všech kolíků.



Chcete-li vložit podložky, postupujte takto:

1. Vyrovnejte podložku s příslušnou pozicí mřížky (A).



- 2. Zatlačte podložku do polohy zastavení.
- Ujistěte se, že podložka se bezpečně zablokovala do segmentu.
 To lze pocítit v posledních několika milimetrech před tím, než se podložka dotkne dorazové polohy.



Chcete-li vyložit podložky, postupujte takto:

- 1. Vytáhněte podložku horizontálně v úrovni plošiny, dokud nebude zcela vyjmuta z plošiny.
- 2. Jednou rukou podepřete přední část podložky.
- 3. Ujistěte se, že se podložka před zvednutím uvolnila ze všech kolíků.



6.3.7 Kontrola uspořádání plošiny

Ujistěte se, že nosiče, laboratorní vybavení a přístroje instalované na plošině odpovídají uspořádání plošiny definovanému pro danou metodu.

OZNÁMENÍ

Poškození zařízení!

Nesprávné umístění segmentů a laboratorního vybavení na pracovním stole může způsobit havárii ramen.

- Vždy se ujistěte, že fyzická konfigurace plošiny a vložené laboratorní vybavení odpovídají konfiguraci pracovního stolu FluentControl.
- Vždy se ujistěte, že je laboratorní vybavení v soupravách správně uloženo. Viz "Pozice laboratorního vybavení" [> 167].

OZNÁMENÍ

Magnetické pole vytváří rušení!

Silné magnetické pole (se severním pólem nahoře) v místě aspirace může rušit snímač přítomnosti špičky a může vést k neočekávaným chybám (např. **ztráta DiTi**).

 Zajistěte, aby v poloze SBS, která sousedí s polohou aspirace, nebyl umístěn žádný silný magnet.



Segmenty Fluent ID může odebrat pouze FSE z důvodu připojení k elektronické desce pod plošinou.

- ✓ Metodu musí připravit hlavní operátor.
- Spotřební materiál odpovídá spotřebnímu materiálu definovanému v metodě.
- Podložky zkumavek Fluent ID musí být vloženy až po spuštění metody na základě výzvy na dotykové obrazovce.
- 1. Postupujte podle pokynů zobrazených na dotykové obrazovce.





Obrázek ukazuje příklad pokynu (A) zobrazeného na dotykové obrazovce:

6.4 Před spuštěním metody

Před spuštěním metody je potřeba vyplnit následující kontrolní seznam.

Přístroj/komponenta	Úkol	
Tab. 15: Kontroly před spuštěním metody		

Přístroj/komponenta	Úkol	Reference/aktivity
Validace procesu	Před zahájením výroby se ujistěte, že zvolená metoda byla validována.	Další informace získáte u hlavního operátora.
Dotyková obrazovka	Postupujte podle pokynů na dotykové obrazovce. OZNÁMENÍ Je potřeba důsledně dodržovat pokyny hlavního operátora.	_
	Pokud se nezobrazí žádné pokyny, postupujte podle níže uvedeného seznamu úloh.	



Přístroj/komponenta	Úkol	Reference/aktivity
Segmenty, nosiče, doplňky a zařízení	Zkontrolujte, zda jsou nainstalovány a zajištěny všechny segmenty, nosiče, doplňky a zařízení.	Pokud se testování nezdaří, obraťte se na hlavního operátora, aby testování provedl znovu.
	Zajistěte, aby se na plošině nacházely pouze předměty určené k použití při této metodě.	
	Zkontrolujte, zda byl test úspěšně dokončen.	
Vzorky a činidla	Zajistěte, aby byly všechny vzorky, činidla a laboratorní vybavení byly správně vloženy.	_
	OZNÁMENÍ Čtení čárových kódů probíhá až po spuštění metody. Před zahájením metody se ujistěte, že na plošině Fluent ID nejsou žádné podložky. Podložky je potřeba vložit, až když se na dotykové obrazovce zobrazí výzva.	
Odpadní potrubí (pouze kapalinové systémy)	Ujistěte se, že odpadní potrubí je vedeno správně.	Vizuálně zkontrolujte, zda není odpadní potrubí zalomené nebo zmačkané. Vyměňte vadné odpadní potrubí. Viz část
		"Připojení mycí stanice (MCA 384)" [▶ 166].
Mycí systém (pouze kapalinové systémy)	Zkontrolujte, zda systémová kapalina a nádoba na odpad jsou správně připojeny.	Viz část "Kontrola potrubí na nádobě systémové kapaliny a nádobě na odpad" [▶ 91].
Mycí systém (pouze kapalinové systémy)	Ujistěte se, že je nádoba systémové kapaliny naplněna na správnou úroveň. Ujistěte se, že je nádoba na odpad prázdná.	Viz část "Připojení nádoby systémové kapaliny a nádoby na odpad" [▶ 142].



Přístroj/komponenta	Úkol	Reference/aktivity
Mycí systém (pouze kapalinové systémy)	Ujistěte se, že je použita správná systémová kapalina, jak je definováno v metodě.	_
Mycí systém (pouze mycí centrum MCA)	Zkontrolujte hladinu kapaliny v mycím bloku.	-
Jednotka odpadu z jednorázových špiček a mycí stanice	Ujistěte se, že jednotka odpadu z jednorázových špiček a mycí stanice je čistá.	Viz část "Čištění jednotky odpadu z jednorázových špiček a mycí stanice" [▶ 136].
	Zajistěte, aby byly namontovány kryty odpadních žlabů pro zadržování aerosolů nebo pro vedení špičky MCA 384.	_
Jednorázové špičky	Zkontrolujte, zda jsou vloženy správné špičky. Ujistěte se, že je odpadní nádoba na špičky prázdná.	_
Fixní hroty	Zkontrolujte, zda jsou fixní hroty čisté a nepoškozené.	Vizuálně zkontrolujte, zda jsou fixní hroty čisté. Vizuálně zkontrolujte fixní hroty pomocí zubního zrcátka, zda je povlak neporušený.
Plošina	Ujistěte se, že nosiče, laboratorní vybavení a přístroje instalované na plošině odpovídají uspořádání plošiny definovanému pro danou metodu.	Viz část "Kontrola uspořádání plošiny" [▶ 87].
Laboratorní vybavení	Zajistěte bezpečné umístění veškerého laboratorního vybavení. Pokud mikrodesky vykazují pohyb do stran, zkontrolujte, zda jsou polohovadla laboratorního vybavení správně nastavená.	Viz část "Pozice laboratorního vybavení" [▶ 167].



Přístroj/komponenta	Úkol	Reference/aktivity
Rotátor zkumavek	Ujistěte se, že na rotátoru zkumavek nechybí žádné polohovací nebo blokovací kolíky.	Viz část "Výměna blokovacích kolíků a polohovacích kolíků" [▶ 194]

6.4.1 Kontrola potrubí na nádobě systémové kapaliny a nádobě na odpad

A UPOZORNĚNÍ

Kontaminace vzorku!

V případě duální kapalinové konfigurace FCA lze pro každé rameno použít různé typy systémových kapalin. Připojení nesprávné nádoby systémové kapaliny k ramenu může způsobit kontaminaci vzorku.

- Označte každou nádobu systémové kapaliny odpovídajícím názvem systémové kapaliny.
- ✓ Mycí systém musí být správně nainstalován.
- 1. Zkontrolujte, zda je hadička (B) správně připojena k nádobě systémové kapaliny (D).
- 2. Pokud je přítomen, zkontrolujte, zda je snímač detekce kapaliny (C) správně připojen k nádobě systémové kapaliny (D).







6.4.2 Kontrola trubek kontejnerů na odpad

1. Zkontrolujte, zda je trubka (A) správně připojena k nádobě na odpad (D).



- 2. Zkontrolujte, zda jsou trubky (C) a (D) správně připojeny k nádobě na odpad.
- 3. Pokud je přítomen, zkontrolujte, zda je snímač detekce kapaliny (E) správně připojen k nádobě na odpad.
- 4. Našroubujte víko (F).



6.5 Spuštění metody

Metoda je soubor skriptů nebo procesů definovaných v softwaru FluentControl. Metodu lze provést v rámci běhu.



Hlavní operátor zapíše metodu, kterou lze provést, následujícím způsobem.

OZNÁMENÍ

Poškození přístroje!

K poškození přístroje může dojít při nesprávném nastavení plošiny nebo při nesprávné obsluze či nesprávném použití softwaru.

- Zkontrolujte, zda jsou všechna bezpečnostní zařízení nainstalována a funkční.
- Ujistěte se, že nosiče, laboratorní vybavení a přístroje instalované na plošině odpovídají uspořádání plošiny definovanému pro danou metodu.
- Zajistěte, aby se na plošině nacházely pouze předměty určené k použití při této metodě.

6.5.1 Spuštění metody

- ✓ Systém FluentControl je spuštěn.
- ✓ Byla provedena část "Přihlášení uživatele" [▶ 81].
- ✓ Byla provedena část "Před spuštěním metody" [▶ 88].
- Vyberte možnost Method Starter (Spuštění metody) (A). Tlačítko se rozsvítí, jakmile se ho dotknete.



Vyberte metodu (B), která se má provést.
 Vybraná metoda je zvýrazněna.



3. Stiskněte tlačítko OK.



4. Stiskněte tlačítko Run (Spustit) (C).

Δ Method Starter	Go 🛞	<
		— C
	Assay 1 is ready to be sta	arted.
		X

- 5. Postupujte podle pokynů na dotykové obrazovce.
- Pokud váš skript obsahuje funkci DeckCheck, věnujte pozornost případným rozdílům ve skutečném živém uspořádání plošiny v porovnání s očekávaným referenčním uspořádání plošiny. Viz "Provoz DeckCheck" [▶ 104].
- 7. Počkejte na ukončení běhu metody.





Na obrazovce se zobrazí přibližný čas (D), kdy bude běh metody ukončen.

8. Pokud se zobrazí hlášení **Run finished with error(s)!** (Cyklus dokončen s chybou!), stiskněte tlačítko **Log** (Protokol) a zkontrolujte chyby a varování.





11:59	inished w	lith er	r or(s)! 12:1
Comment 2 E: Script execution t 1 Run paused.	Process erminated by End Script star	Log	Evd Ter 12:00:48 (2017-08-18) 11:59:35 (2017-08-18)

9. Stisknutím tlačítka Next (Další) se vrátíte na úvodní obrazovku.

10. Stiskněte tlačítko **Exit** (Konec).

Po ukončení běhu metody se zobrazí následující obrazovka (E).

۵	Running 🛞	
		_
	Run finished successfully.	
	10:09 10: Start Time End 1	43 Ime
L.		_
	E	
		1
Show		Exit



D UPOZORNĚNÍ



Produkt Fluent je laserový výrobek třídy 1 podle normy IEC 60825-1:2014, který vyzařuje laserové záření.

Laserový paprsek může způsobit oslnění, zábleskovou slepotu a paobrazy.Nedívejte se do laserového paprsku ani do jeho odrazů.



6.5.2.1 Vkládání podložek Fluent ID

OZNÁMENÍ

Poškození v důsledku nesprávného vkládání nebo vykládání

Poškození podložek a kolíků.

- Vyrovnejte podložku horizontálně s plošinou.
- Jednou rukou podepřete přední část podložky.
- Při vykládání se ujistěte, že se podložka před zvednutím uvolnila ze všech kolíků.
- ✓ Produkt Fluent je vybaven skenerem čárových kódů Fluent ID.
- ✓ Zkumavky se vkládají do podložek štítkem s čárovým kódem směrem doleva.
- ✓ Všechny zkumavky v podložce mají stejnou velikost a tvar. Typy podložek zkumavek uvádí část "Podložky pro kompatibilní zkumavky Fluent ID" [▶ 60].
- 1. Zvolte a spusťte metodu pomocí dotykové obrazovky.

LED kontrolky začnou blikat a objeví se zpráva **Prosím, vložte zkumavky** na dotykové obrazovce.

Při použití různých typů zkumavek dbejte na to, aby byl pro každou konkrétní mřížku použit správný typ podložky.

- 2. Jednou rukou podepřete přední část podložky.
- 3. Podržte podložku ve vodorovné poloze na úrovni plošiny.
- 4. Zatlačte podložku do polohy zastavení.







5. Posuňte podložky zkumavek jednu po druhé na vyhrazené mřížky v oblasti pro vkládání Fluent ID.

Zkontrolujte, zda byly úspěšně naskenovány všechny čárové kódy.
 LED kontrolky se rozsvítí zeleně, když jsou podložky zkumavek ve vložené poloze a všechny štítky s čárovým kódem byly úspěšně naskenovány.

Popis stavu LED Fluent ID naleznete v části "Stavové LED kontrolky Fluent ID" [> 77].



- 7. V případě chyby při snímání čárového kódu vyjměte podložku, napravte problém a znovu podložku vložte.
- 8. Táhněte podložku vodorovně podél plošiny, dokud není zcela vyjmuta.





6

Fluent ID přečte každý kód při průchodu skenerem několikrát. U malých a úzkých zkumavek (tj. průměr ≤ 10 mm) snižte rychlost ručního vkládání, abyste umožnili všechna čtení a snížili počet chybových hlášení.



Errors: 🔲 Missing Runner 🗖 Wrong Runner 🖲 Unreadable 🍨 Duplicates O Missing

Obr. 42: Potvrzení načtení čárového kódu zobrazené na dotykovém displeji

Tab. 16: Význam GUI (podložka)

Obdélník (podložka)	Význam
Zelená	Všechny čárové kódy zkumavek v podložce byly úspěšně načteny.
Bílá s červeným okrajem	Chybný typ podložky pro tuto pozici na mřížce.
Šedá s červeným okrajem	Chybějící podložka. Do této pozice mřížky by měla být vložena podložka.

Tab. 17: Význam GUI (pozice zkumavky)

Kroužek (pozice zkumavky)	Význam
Zelená	Čárové kódy byly úspěšně načteny.
Červená	Nečitelný čárový kód
Oranžová	Duplicitní čárový kód
Bílá s červeným okrajem	Chybějící zkumavka. Do této pozice by měla být vložena zkumavka.

6

Při použití podložky pro 2 ml zkumavky Safe-Lock není možné odlišit chybějící zkumavky od nečitelných čárových kódů. Chybějící zkumavky jsou hlášeny jako nečitelné čárové kódy.

6.5.2.2 Vykládání podložek Fluent ID

- Běh byl dokončen nebo běh právě probíhá a kontrolky LED blikají se zprávou
 Prosím, vyložte zkumavky zobrazenou na dotykové obrazovce.
- 1. Táhněte podložku vodorovně podél plošiny, dokud není zcela vyjmuta.



6.5.3 Vkládání a vykládání podložek rotátoru zkumavek

6.5.3.1 Vkládání podložek rotátoru zkumavek

DEPOZORNĚNÍ

Biologická kontaminace systému a/nebo uživatele!

Poškozené zkumavky na vzorky mohou implodovat, což může vést k rozlití vzorku na rotátor zkumavek.

- Dbejte, aby do rotátoru zkumavek nebyly vkládány poškozené zkumavky.
- ✓ Produkt Fluent je vybaven rotátorem zkumavek.
- Zkumavky se vkládají do podložek rotátoru zkumavek se štítkem s čárovým kódem směrem doleva.
- ✓ Všechny zkumavky v podložce mají stejnou velikost a tvar. Typy podložek zkumavek uvádí část "Podložky rotátoru zkumavek" [▶ 63].
- 1. Zvolte a spusťte metodu pomocí dotykové obrazovky.

LED kontrolky začnou blikat a objeví se zpráva **Prosím, vložte zkumavky** na dotykové obrazovce. Při použití různých typů zkumavek se ujistěte, že jste zvolili správnou podložku pro daný typ zkumavky (BD nebo Greiner). Dbejte také na to, abyste do příslušných podložek vkládali zkumavky různých výšek: Zkumavky jsou vždy drženy ve své poloze podložkou ve výšce jejich uzávěrů. Dna zkumavek musí vždy pevně sedět ve vložkách zkumavek na podložkách.

2. Otevřete zajišťovací páku podložky.







3. Jednou rukou podepřete přední část podložky.

- 4. Podržte podložku ve vodorovné poloze na úrovni plošiny.
- 5. Zatlačte podložku do polohy zastavení.
- 6. Na vyhrazené mřížky na rotátoru zkumavek postupně nasuňte podložky rotátoru zkumavek.

$ \Delta angle$ Running $ angle$ Load tube	runr	ners	ŧ.					
	000000000000000000000000000000000000000	000000000000000000000000000000000000000	000000000000000000000000000000000000000	000000000000000000000000000000000000000	000000000000000000000000000000000000000	000000000000000000000000000000000000000		
	1	2	3	4	5	6		
Please load tube runners a	nd pres	is con	tinue v	when	finish	ned	Errors: 🔲 Mi	ssing Runner
							-	~
							Stop	Continue

- Zkontrolujte, zda byly úspěšně naskenovány všechny čárové kódy. Kontrolky LED se rozsvítí zeleně, když jsou podložky zkumavek ve vložené poloze a všechny štítky s čárovým kódem byly úspěšně naskenovány. Popis stavu LED diody rotátoru zkumavek naleznete v části "Stavové LED kontrolky Fluent ID" [▶ 77].
- 8. V případě chyby při snímání čárového kódu vyjměte podložku, napravte problém a znovu podložku vložte.





9. Zavřete zajišťovací páku podložky.

00000000000		000000000000000000000000000000000000000	000000000000000000000000000000000000000	000000000000000000000000000000000000000	000000000000
1	2	3	4	5	6

Errors: Missing Runner Wrong Runner • Unreadable • Duplicates • Missing Obr. 43: Potvrzení načtení čárového kódu zobrazené na dotykovém displeji

Tab. 18: Význam GUI (podložka)

Obdélník (podložka)	Význam
Zelená	Všechny čárové kódy zkumavek v podložce byly úspěšně načteny.
Bílá s červeným okrajem	Chybný typ podložky pro tuto pozici na mřížce.
Šedá s červeným okrajem	Chybějící podložka. Do této pozice mřížky by měla být vložena podložka.

Tab. 19: Význam GUI (pozice zkumavky)

Kroužek (pozice zkumavky)	Význam		
Zelená	Čárové kódy byly úspěšně načteny.		



Kroužek (pozice zkumavky)	Význam
Červená	Nečitelný čárový kód
Oranžová	Duplicitní čárový kód
Bílá s červeným okrajem	Chybějící zkumavka. Do této pozice by měla být vložena zkumavka.

6.5.3.2 Vykládání podložek rotátoru zkumavek



Neukládejte podložky rotátoru zkumavek s vloženými zkumavkami jinak než v provozních podmínkách manipulace s kapalinami. Viz část "Podmínky prostředí" [> 44].

- Běh byl dokončen nebo běh právě probíhá a kontrolky LED blikají se zprávou Prosím, vyložte zkumavky zobrazenou na dotykové obrazovce.
- Rotátor trubek je v horizontální domovské poloze.
- 1. Otevřete zajišťovací páku podložky.



2. Táhněte podložku vodorovně podél plošiny, dokud není zcela vyjmuta.



Segmenty rotátoru zkumavek může odebrat pouze FSE z důvodu připojení k elektronické desce pod plošinou.

- Metodu musí připravit hlavní operátor.
- Spotřební materiál odpovídá spotřebnímu materiálu definovanému v metodě.
- Podložky rotátoru zkumavek musí být vloženy až po spuštění metody na základě výzvy na dotykové obrazovce.



6.5.4 Resetování chyb

Pokud se zobrazí hlášení, postupujte následovně:

Hlášení

- Zkontrolujte funkci displeje, funkci tlačítek nebo chybové hlášení. Viz části "Pracovní oblast" [▶ 71] a "Tlačítka obnovení metody" [▶ 72].
- 2. Při opravě chyby postupujte podle pokynů v této příručce a na dotykovém displeji.
- Pokračujte v běhu metody. Viz část "Tlačítka zobrazení, možností a akcí" [▶ 71].

Pokud se stavová kontrolka rozsvítí nebo změní barvu, postupujte následovně:

Stavová kontrolka

- 1. Zkontrolujte stav přístroje. Viz část.
- Je-li produkt Fluent vybaven snímačem čárových kódů Fluent ID, zkontrolujte stav LED snímače čárových kódů Fluent ID. Viz část "Stavové LED kontrolky Fluent ID" [▶ 77].
- 3. Zkontrolujte funkci displeje, funkci tlačítek nebo chybové hlášení. Viz části "Pracovní oblast" [▶ 71] a "Tlačítka obnovení metody" [▶ 72].
- 4. Podívejte se do tabulky řešení problémů. Viz část "Tabulky pro řešení problémů" [▶ 152].
- 5. Pokud problém nelze vyřešit, viz "Zákaznická podpora" [▶ 199].

6.6 Provoz DeckCheck

Pokud váš skript zahrnuje použití systému DeckCheck, kamerový systém DeckCheck po načtení vyfotí pracovní stůl a porovná skutečné živé uspořádání s referenčním uspořádáním.

DeckCheck potřebuje přibližně 20 sekund v případě systému se třemi rameny / třemi kamerami a přibližně 12 sekund pro systém s jedním nebo dvěma rameny / jednou kamerou, aby pořídil snímky plošiny a zobrazil srovnání živého a referenčního uspořádání (za předpokladu, že konfigurace počítače je správná – viz příručku aplikačního softwaru FluentControl).

Všimněte si, že pouze při prvním použití po zapnutí přístroje bude příkaz DeckCheck potřebovat delší dobu k zobrazení prvního výsledku – může to trvat několik minut.

Během této doby bude svítit zadní LED dioda.

U tříramenných systémů Fluent se musí prostřední rameno pohybovat mezi levou a pravou polohou (u jednoramenných nebo dvouramenných systémů bude levé a pravé rameno umístěno na krajní levé, resp. krajní pravé straně). Pro tento pohyb ramena musí být přední dveře zavřené. Pokud je snímek pořizován při otevřených dveřích, jedna kamera je zpravidla blokována prostředním ramenem nebo jiným ramenem, které bylo ručně posunuto.

Během procesu DeckCheck se na dotykové obrazovce zobrazují tvary, které se pohybují na obrazovce a dále **Taking Images** po kterém následuje **Checking**. Po 12-20 sekundách se v závislosti na velikosti a konfiguraci přístroje zobrazí obrázky plošiny ve střídavém režimu. Zobrazí se obraz první kamery se zaznamenanou odchylkou počínaje zleva.

Obrazovka DeckCheck zobrazující nesrovnalost ve vztahu k referenčnímu obrázku. Zde je na referenčním snímku vidět, že by měla být přítomna plošina a středová kamera zjistila nesrovnalost.





Pokud je vám nabídnuta možnost **Ignore & Continue** dole, běh bude pokračovat s dominantním pracovním stolem. Stiskněte tlačítko **Ignore & Continue** v případě, že jste si jisti, že již nejsou žádné rozdíly oproti požadovanému pracovnímu stolu a před zavřením dvířek. Vyberte **Check**, pokud chcete pořídit nové snímky plošiny – mějte na paměti, že pokud nejsou dveře u tříramenných systémů zavřené, obraz bude pořizován, ale prostřední rameno bude kameře překážet. Kontrola se jinak provádí automaticky při zavření dveří.



Všimněte si, že pokud skript obsahuje možnost **vždy zobrazit**, zobrazí se výše uvedená obrazovka a nejsou zvýrazněny žádné nesrovnalosti. Referenční a živý obraz se však budou střídat a může se stát, že se objeví malé rozdíly, které systém nezachytí, ale jsou snadno viditelné okem – například některé barevné rozdíly, chybějící zkumavky nebo malé boční posuny. Viz limity uvedené níže.

Pokud jsou zjištěny nesrovnalosti, budou zvýrazněny.

Chcete-li opravit rozdíly:



- 1. Otevřete dvířka a vyměňte nebo opravte polohu zvýrazněných položek.
- 2. DeckCheck bude průběžně porovnávat opravenou živou situaci s referenčním uspořádáním.
- Pomocí tlačítek DeckCheck se můžete podívat na rozdíly zachycené jednotlivými kamerami nebo podle potřeby pozastavit zobrazení a podržet referenční obraz. Pokud nejsou zjištěny žádné další rozdíly, zobrazí se zelené tlačítko pokračovat.
- 4. Vyberte Continue pro pokračování v metodě.



Pokud jsou zbývající rozdíly považovány za přijatelné (např. celkový počet špiček může být na začátku metody proměnlivý nebo hladiny kapaliny se na začátku běhu výrazně liší), můžete zvolit možnost **Ignore & Continue** v případě, že ji hlavní operátor ve skriptu nabídne.

Na některé rozdíly v uspořádání nemusí DeckCheck upozornit – např. na následující barevné zásobníky špiček FCA:

Rozdíl mezi typy adaptérů hlavy MCA:

- Žlutá/oranžová
- Bílá/oranžová
- Šedá / všechny barvy

MCA 384, různé typy špiček

MCA 96, různé typy špiček

Chybějící zkumavky na částečně naplněných podložkách zkumavek

Žlábek 300 SBS

Mikrodesky otočené o 180 stupňů

Tvar jamky mikrodesky (např. kulatá nebo s plochým dnem nebo jamka PCR)

Desky v periferních hotelech 10 ml/ 25 ml, žlábky jako vložka

Částečně průhledná víka

Mnohé z těchto rozdílů jsou však jasně patrné při přepínání mezi živým a referenčním uspořádáním.

6.7 Obnovení metody

FluentControl nabízí možnost obnovení po chybách - např.:

Předchozí běh metody byl přerušen nebo došlo k fatální chybě: Možnost obnovení metody nabízí možnost pokračovat od bodu, ve kterém došlo k selhání v předchozím běhu.



Po přerušení metody nebo fatální chybě musí být provedena denní údržba. Viz "Denní péče o systém" [> 114].



6.7.1 Přepnutí do režimu obnovení metody

- ✓ Hlavní operátor povolil ve FluentControl možnost obnovení metody.
- ✓ Předchozí běh metody byl přerušen.
- 1. Vyberte Obnovení metody (A).



6.7.2 Obnovení běhu metody

- ✓ Část "Přepnutí do režimu obnovení metody" [▶ 107] byla provedena.
- 1. Vyberte možnost pokračování na další obrazovku (A).

Na obrazovce se zobrazí poslední provedený řádek skriptu (C) a řádek skriptu, kde došlo k selhání – bod obnovení (B).

Δ Method Recov	ery 🛞
E Reagent Distribu	tion
Sample Transfer	— C
Get Tips	PCA 1 PCA, 200ul 585
Aspirate	FCA 1 100 (µl) Water Video Single 1x24 13x75mm Tube Runner(001) Q1 - X1
Uispense	FCA t 100 [µl] Water Video Single 96 Well Flat[001] A1 - H1
Drop Tips	FCA 1 Wash Station Thru Deck DiTi Waste_1
🗐 🏣 Incubate & Wash	
Read & Export D	ata
•	A → ×
Discard	Worktable Cancel

 Vyberte požadovaná tlačítka popsaná v části "Tlačítka obnovení metody" [▶ 72].



- 3. Ujistěte se, že fyzické rozložení plošiny produktu Fluent odpovídá rozložení pracovního stolu (D) zobrazenému na dotykové obrazovce.
- 4. Vyberte Obnovení běhu (E).

Systém se spustí.



6.8 Vypnutí přístroje

Pokud není spuštěna žádná metoda, přístroj se přepne do pohotovostního režimu. Přístroj v elektrické síti není nutné vypínat.

Chcete-li přístroj vypnout, postupujte následovně:

- 1. Zastavte metodu a na dotykovém displeji zvolte pohotovostní režim.
- 2. Zkontrolujte, zda je přístroj v pohotovostním režimu. Viz část .
- 3. Umístěte robotická ramena do volného prostoru pr pohyb.






4. Vypněte napájení vypínačem (A) na zadní straně externího zdroje napájení.



7 Péče o systém

V této kapitole jsou uvedeny pokyny pro všechny úkoly péče o systém, které je třeba provést pro udržení přístroje Fluent v dobrém provozním stavu.



Přístroj Fluent používejte, pouze pokud je v dobrém provozním stavu. Důsledně dodržujte pokyny pro péči o systém popsané v tomto návodu. Pro zajištění optimálního výkonu a spolehlivosti provádějte pravidelně údržbu a čištění.

V případě jakýchkoli problémů a dotazů se podívejte do části "Zákaznická podpora" [▶ 199].



DEPOZORNĚNÍ

Rameno naráží do objektů na pracovním stole

S rameny Fluent lze pohybovat manuálně. Manuální pohyb ramen musí být prováděn jemně, rameno je nutné držet. Nenarážejte rameny do pevných předmětů, a to ani do mechanické zarážky rozsahu ramen

7.1 Dekontaminace

Dekontaminace se podle standardních laboratorních předpisů vyžaduje za okolností uvedených v části "Prohlášení o dekontaminaci" [> 39].

A VAROVÁNÍ

Kontaminace!

Zbytky látek na přístroji Fluent mohou způsobit úraz a ovlivnit integritu procesu.

 Před každou interakcí proveďte dekontaminaci produktu Fluent a všech jeho dílů a příslušenství.

Metodu dekontaminace musí určit hlavní operátor na základě typu kontaminantu a stupně kontaminace. V této kapitole jsou uvedeny pokyny pro výběr dekontaminačních prostředků a způsobů jejich použití.



Informace o ošetření parami peroxidu vodíku naleznete v Referenční příručce. Viz Referenční dokumenty.

A UPOZORNĚNÍ

Nesprávné výsledky měření pomocí čtečky Frida!

Pokud není namontována žádná vložka, může čtečka Frida poskytovat nesprávné výsledky měření.

 Pokud je vložka vyjmuta (např. kvůli čištění), použijte červenou zaslepovací zátku.



7.2 Čisticí prostředky

7.2.1 Specifikace čisticích prostředků

Pro péči o systém jsou nutné speciální čisticí prostředky. Všechny doporučené čisticí prostředky byly pečlivě vybrány a testovány.

OZNÁMENÍ

Snížená účinnost a chemická kompatibilita!

Účinnost čisticích prostředků a chemická kompatibilita nejsou zaručeny, pokud jsou použity jiné čisticí prostředky než ty, které doporučuje společnost Tecan.

- Používejte pouze čisticí prostředky doporučené společností Tecan.
- Čisticí prostředky jsou pro každé konkrétní použití definovány v tabulkách péče o systém. Nepoužívejte čisticí prostředky, pokud nejsou určeny k použití pro konkrétní úkol.

Následující tabulka specifikuje čisticí prostředky určené k použití podle tabulek péče o systém a činností péče o systém:

Činidlo	Specifikace
DI voda	Destilovaná nebo deionizovaná voda
Alkohol	70% etanol, 100% isopropanol (2-propanol)
Slabý detergent	Liqui-Nox
Povrchově aktivní látka	Contrad 70, Contrad 90 / Contrad 2000, Decon 90
Dezinfekční prostředek	Bacillol plus, SporGon
Dezinfekční prostředek na povrchy (na kontaminaci nukleovými kyselinami)	DNAzap
Slabá kyselina	kyselina sírová 0,3M, 10% kyselina octová, 30–40% kyselina mravenčí
Zásada	hydroxid sodný 0,1M
Bělidlo	2% chlornan sodný
Systémová kapalina	Jak je definováno v metodě. Vezměte na vědomí, že vodné roztoky s obsahem solí by se měly během nečinnosti systému – např. přes noc nebo o víkendech – vyplachovat. Viz část Péče o systém – "Konec dne" [▶ 116].

Tab. 20: Čisticí prostředky určené k použití podle tabulek péče o systém



7.2.2 Komerční čisticí prostředky

Je potřeba si pečlivě přečíst a dodržovat všechny pokyny výrobce čisticích prostředků nebo pokyny uvedené v tomto návodu pro manipulaci s čisticími prostředky.

V následující tabulce je uveden seznam komerčně dostupných čisticích a dezinfekčních prostředků určených k použití, jak je uvedeno v tabulkách péče o systém a v činnostech péče o systém.

Čisticí prostředek	Kategorie činidla	Výrobce
DNAzap	Dezinfekční prostředek na povrchy (pro povrchy kontaminované nukleovými kyselinami)	Ambion www.ambion.com
Decon, Contrad	Povrchově aktivní látka	Decon Laboratories www.deconlabs.com
SporGon	Dezinfekční prostředek	Decon Laboratories www.deconlabs.com
Bacillol Plus	Dezinfekční prostředek	www.bode-chemie.com
Liqui-Nox	Slabý detergent	Alconox www.alconox.com

Tab. 21: Komerční čisticí prostředky

7.3 Režim péče o systém

Hlavní operátor definuje požadované metody péče o systém podle tabulek péče o systém uvedených v dokumentu "Tabulky péče o systém" [> 114]. Péče o systém přístupná na dotykové obrazovce uvádí pokyny pro úkoly péče o systém.

7.3.1 Přepnutí do režimu péče o systém

✓ Musí být k dispozici metody péče o systém.



1. Vyberte Péče o systém (A).



2. Vyberte úkol, který se má provést.

☆ System Care ⊗	<
Start of Day End of Day Weekly	
Der is not exceeded Der is not exceeded + +	
T T	
Remove	Cancel

- 3. Stiskněte **Běh** pro iniciaci metody péče o systém.
- 4. Proveďte úkoly péče o systém.

7.3.2 Resetování chyb

Pokud se zobrazí hlášení, postupujte následovně:

Hlášení

- Zkontrolujte funkci displeje, funkci tlačítek nebo chybové hlášení. Viz části "Pracovní oblast" [▶ 71] a "Tlačítka obnovení metody" [▶ 72].
- 2. Při opravě chyby postupujte podle pokynů v této příručce a na dotykovém displeji.
- Pokračujte v běhu metody. Viz část "Tlačítka zobrazení, možností a akcí" [▶ 71].

Pokud se stavová kontrolka rozsvítí nebo změní barvu, postupujte následovně:



Stavová kontrolka

- 1. Zkontrolujte stav přístroje. Viz část.
- Je-li produkt Fluent vybaven snímačem čárových kódů Fluent ID, zkontrolujte stav LED snímače čárových kódů Fluent ID. Viz část "Stavové LED kontrolky Fluent ID" [▶ 77].
- Zkontrolujte funkci displeje, funkci tlačítek nebo chybové hlášení. Viz části "Pracovní oblast" [▶ 71] a "Tlačítka obnovení metody" [▶ 72].
- Podívejte se do tabulky řešení problémů. Viz část "Tabulky pro řešení problémů" [▶ 152].
- 5. Pokud problém nelze vyřešit, viz "Zákaznická podpora" [▶ 199].

7.4 Tabulky péče o systém

Pro zajištění optimálního výkonu a spolehlivosti provádějte pravidelně údržbu a čištění dle doporučení.

6

Úkoly v tabulkách péče o systém lze provádět pouze v režimu péče o systém. Viz část "Režim péče o systém" [▶ 112].

Úkoly péče o systém je třeba provádět v pravidelných intervalech – denně, týdně a měsíčně.

7.4.1 Denní péče o systém

7.4.1.1 Začátek dne

Spusťte **DailySystemCare**, pokud ji hlavní operátor zpřístupní; nebo proveďte jednotlivé úlohy, které se vztahují k vaší konfiguraci ramena Fluent a jsou uvedeny v následující tabulce v chronologickém pořadí.

Tab. 22: Tabulka péče o systém na začátku dne

Přístroj / Komponenta	Úloha péče o systém	Čisticí prostředek / Jednorázový výrobek / Zařízení	Reference / Činnosti péče o systém
Propichovací špičky	Vizuálně zkontrolujte, zda na propichovacích špičkách nejsou usazeniny. V případě potřeby vyčistěte. Zkontrolujte, zda nejsou špičky ohnuté.	70% ethanol nebo 2% bělidlo a hadřík, který nepouští vlákna.	Viz část "Čištění propichovacích špiček" [▶ 127].
Jednorázové kužely špiček a pevné špičky	Kontrola poškození a usazenin		Tento úkol je zahrnut v metodě Denní péče o systém . OZNÁMENÍ Jednorázové špičky nejsou určeny k opakovanému použití.



Přístroj / Komponenta	Úloha péče o systém	Čisticí prostředek / Jednorázový výrobek / Zařízení	Reference / Činnosti péče o systém
Pevné špičky	Vyčistěte. Zkontrolujte, zda nejsou špičky ohnuté. Vizuálně zkontrolujte pomocí zubního zrcátka, zda je povlak neporušený.	70% ethanol nebo 100% isopropanol a hadřík, který nepouští vlákna.	Viz část "Čištění fixních hrotů" [▶ 127].
Nádoba na systémovou kapalinu (Rameno Liquid FCA a MCA s pevnými špičkami)	Ujistěte se, že je čisté a plné bez viditelných bublin Zkontrolujte, zda jsou hadice ke konektorům nádoby správně připojeny.	_	Tento úkol je zahrnut v metodě Denní péče o systém .
Nádoba na tekutý odpad (Rameno Liquid FCA a MCA s pevnými špičkami)	Ujistěte se, že je prázdný Zkontrolujte, zda jsou hadice ke konektorům nádoby správně připojeny.	_	Tento úkol je zahrnut v metodě Denní péče o systém .
Sáček na odpad jednorázových špiček	Ujistěte se, že je prázdný	_	Viz část "Výměna sáčku na odpad jednorázových hrotů" [▶ 139]. Tento úkol je zahrnut v metodě Denní péče o systém .
Kapalinový systém (Liquid FCA)	Ujistěte se, že je čistý	Systémová kapalina, alkohol, DI voda	Tento úkol je zahrnut v metodě Denní péče o systém nebo může být spuštěn samostatně jako metoda Rutinní údržba proplachování ramena Liquid FCA . Viz část "Čištění cesty kapaliny" [▶ 141].
Mycí systém (MCA s pevnými špičkami)	Proplach/plnění	Mycí kapalina	Spusťte přímý příkaz Plnicí mycí stanice (MCA384) .
Kapalinový systém (Liquid FCA)	Po propláchnutí vizuálně zkontrolujte, zda na špičkách nebo kuželu DiTi nejsou kapky	_	Tento úkol je zahrnut v metodě Denní péče o systém .



Přístroj / Komponenta	Úloha péče o systém	Čisticí prostředek / Jednorázový výrobek / Zařízení	Reference / Činnosti péče o systém
Prsty upínadla	Zkontrolujte, zda jsou prsty přímé a vyrovnané Kontrola poškození a chybného vyrovnání		V případě chybného vyrovnání viz část "Řešení problémů s robotickým upínacím ramenem (RGA – Robotic Gripper Arm)" [▶ 162]. Deformace nebo poškození. Viz část "Zákaznická podpora" [▶ 199].
Prsty upínadla FCA	Kontrola poškození	_	V případě poškození proveďte výměnu. Informace o objednávání naleznete v Referenční příručce. Viz část Referenční dokumenty.
Čtečka Frida	Vyjměte zaslepovací zátku a nasaďte vložku do čtečky Frida	_	Viz část "Čtečka Frida" [▶ 149].

A UPOZORNĚNÍ

Nesprávné výsledky měření pomocí čtečky Frida!

Pokud není namontována žádná vložka, může čtečka Frida poskytovat nesprávné výsledky měření.

 Pokud je vložka vyjmuta (např. kvůli čištění), použijte červenou zaslepovací zátku.

7.4.1.2 Konec dne

V následující tabulce jsou uvedeny úkoly denní péče o systém na konci dne v chronologickém pořadí:

Tab.	23:	Tabulka	péče c	systém	na	konci	dne
------	-----	---------	--------	--------	----	-------	-----

Přístroj / Komponenta	Úloha péče o systém	Čisticí prostředek / Jednorázový výrobek / Zařízení	Reference / Činnosti péče o systém
Podnosy pracovní plochy	Zkontrolujte rozlité kapaliny a podle potřeby je vyčistěte nebo vyměňte.	DI voda, alkohol, slabý čisticí prostředek, dezinfekční prostředek, zásada, bělidlo, DNAzap	Viz část "Čištění podnosů plošiny" [▶ 132].



Přístroj / Komponenta	Úloha péče o systém	Čisticí prostředek / Jednorázový výrobek / Zařízení	Reference / Činnosti péče o systém
Segmenty Kryt Fluent ID	Vyčistěte	DI voda, alkohol, slabý čisticí prostředek, dezinfekční prostředek, zásada, bělidlo, DNAzap OZNÁMENÍ Okno skeneru vyžaduje jiné čisticí prostředky než	Viz část "Čištění podložek a segmentů" [▶ 132]. VAROVÁNÍ! Nedívejte se do laserového paprsku.
		samotný segment. Viz "Týdenní péče o systém" [▶ 119].	
Reflexní fólie (Fluent ID, rotátor zkumavek)	Vyčistěte a zkontrolujte, zda nedošlo k poškození	Alkohol OZNÁMENÍ Reflexní fólie vyžaduje jiné čisticí prostředky než samotný segment.	Poškození. Viz část "Výměna reflexní fólie Fluent ID" [▶ 134].
Podložky	Vyčistěte	DI voda, alkohol, slabý čisticí prostředek, dezinfekční prostředek, povrchově aktivní látka, slabá kyselina, zásada, bělidlo, DNAzap	Viz část "Čištění podložek a segmentů" [▶ 132].
Blok pevných špiček (MCA)	Zaparkujte, vyčistěte, zkontrolujte a zakryjte	Alkohol	Viz část "Čištění bloku fixních špiček (MCA 384)" [▶ 147]. UPOZORNĚNÍ! Společnost Tecan doporučuje na konci každé pracovní směny blok špiček zaparkovat a vyjmout z přístroje. Blok špiček vždy vyčistěte a náležitě uložte. Zkontrolujte posunutí špičky.
Pevné špičky	Vyčistěte	Alkohol, bělidlo, hadřík bez žmolků	Viz část "Čištění fixních hrotů" [▶ 127].
Kužel na jednorázové špičky	Vyčistěte	Alkohol, hadřík bez žmolků	Viz část "Čištění jednorázových kuželů hrotů" [▶ 126].



Přístroj / Komponenta	Úloha péče o systém	Čisticí prostředek / Jednorázový výrobek / Zařízení	Reference / Činnosti péče o systém
Mycí a odpadní stanice (Liquid FCA)	Vyčistěte	DI voda, alkohol, slabý čisticí prostředek, dezinfekční prostředek	Viz část "Čištění jednotky odpadu z jednorázových špiček a mycí stanice" [▶ 136].
Skluz odpadních jednorázových špiček a odpadní kryty	Vyčistěte	DI voda, alkohol, slabý čisticí prostředek, dezinfekční prostředek	Viz část "Čištění skluzu odpadních jednorázových hrotů" [▶ 137].
Kapalinový systém (Liquid FCA)	Propláchněte	Systémová kapalina OZNÁMENÍ Pokud má kapalinový systém vysoký obsah soli, propláchněte jej deionizovanou vodou.	Spusťte metodu Liquid FCA Routine Flush Maintenance (Údržba proplachování ramena Liquid FCA).
Mycí stanice MCA (MCA s pevnými špičkami)	Vyčistěte	Alkohol, bělidlo, Decon 90, Contrad OZNÁMENÍ Proces mohou ovlivnit povrchově aktivní látky, jako je Decon nebo Contrad. Pokud tedy takové prostředky používáte, pečlivě proces validujte.	Spusťte přímý příkaz Prime Wash Station (MCA384).
Sáček na odpad jednorázových špiček	Vyměňte	Doporučené specifikace sáčku: Š × D: 300 mm × 600 mm; tloušťka: 0,5 mm Materiál: Polypropylen, polyethylen nebo kopolymer (s možností ošetření v autoklávu) OZNÁMENÍ Použitý sáček na odpad musí být v souladu s místními bezpečnostními předpisy.	Viz část "Čištění skluzu odpadních jednorázových hrotů" [▶ 137].



Přístroj / Komponenta	Úloha péče o systém	Čisticí prostředek / Jednorázový výrobek / Zařízení	Reference / Činnosti péče o systém
Nádoba na systémovou kapalinu (Liquid FCA)	Ujistěte se, že je čistý	Systémová kapalina	Viz část "Připojení nádoby systémové kapaliny a nádoby na odpad" [▶ 142].
Odpadní nádoba (Rameno Liquid FCA a MCA s pevnými špičkami)	Vyprázdnění a vyčištění	Deionizovaná voda, alkohol, slabý čisticí prostředek, povrchově aktivní látka, dezinfekční prostředek, zásada, bělidlo	V závislosti na místních laboratorních předpisech čistěte denně nebo týdně. Viz část "Připojení nádoby systémové kapaliny a nádoby na odpad" [▶ 142].
Bezpečnostní panel	Vyčistěte	DI voda, alkohol, slabý čisticí prostředek	Viz část "Čištění bezpečnostních panelů" [▶ 136].
Rotátor zkumavek	Vyčistěte povrchy, držák a mycí stanici	Hadříky bez žmolků s 2% bělidlem, 70% etanolem nebo 100% isopropanolem.	Viz část "Čištění rotátoru zkumavek" [▶ 127].
Propichovací špičky	Vizuálně zkontrolujte, zda na propichovacích špičkách nejsou usazeniny. V případě potřeby vyčistěte.	70% ethanol nebo 2% bělidlo a hadřík, který nepouští vlákna.	Viz část "Čištění propichovacích hrotů" [▶ 127].
Mycí stanice – propichovací špičky a odpadní hadičky	Vyčistěte mycí stanici a hadičky od zbytků vzorku.	DI voda, slabý čisticí prostředek, dezinfekční prostředek. Nepoužívejte bělidlo, pokud následně nepropláchnete součásti mycí stanice vodou.	Proveďte prodloužený proplach mycí stanice, a to všech jejích částí (přední, střední i zadní části).
Čtečka Frida	Vyjměte vložku a nasaďte zaslepovací zátku do čtečky Frida.	-	Viz část "Čtečka Frida" [▶ 149].

7.4.2 Týdenní péče o systém

Týdenní péče o systém by se měla provádět v poslední pracovní den každého týdne.

Spusťte **metodu WeeklySystemCare**, pokud ji hlavní operátor zpřístupní; nebo navíc k denním úkolům proveďte jednotlivé úlohy, které se vztahují k vaší konfiguraci ramena Fluent a jsou uvedeny v následující tabulce v chronologickém pořadí.



Tab. 24: Týdenní tabulka péče o systém

Přístroj / Komponenta	Úloha péče o systém	Čisticí prostředek / Jednorázový výrobek / Zařízení	Reference / Činnosti péče o systém
Mycí systém (MCA 384 s pevnými špičkami)	Vizuálně zkontrolujte, zda nejsou v hadicích a filtru nečistoty	Filtr	Znečištěný filtr. Viz část "Zákaznická podpora" [▶ 199].
Pipetovací hlava (MCA 384)	Proveďte zkoušku těsnosti	_	Spusťte metodu kontroly těsnosti a nulového dávkování ramena MCA 384
Adaptér desky (MCA 384)	Vyčistěte	Alkohol Stlačený vzduch (pro sušení)	Viz část "Čištění adaptéru desky (MCA 384)" [▶ 147].
Kapalinový systém (Liquid FCA)	Vyčistěte	V závislosti na kapalině, kterou zpracovává Fluent Decon, Contrad, zásada, slabá kyselina, dezinfekční prostředek Následují proplachy vodou, alkoholem a systémovou kapalinou	Viz část "Čištění cesty kapaliny" [▶ 141].
Liquid FCA	Zkontrolujte správnou těsnost stříkaček na rozhraní ventilu a správnou těsnost pístu stříkačky na zajišťovacím šroubu pístu.	_	Viz část "Kontrola těsnosti stříkaček" [▶ 142]
Kužely DiTi	Kontrola těsnosti kuželu DiTi	_	Viz část "Utahování kužele DiTi" [▶ 148]
Liquid FCA	Proveďte zkoušku těsnosti (Liquid FCA)	-	Spusťte zkoušku těsnosti ramena Liquid FCA.
Air FCA	Proveďte zkoušku těsnosti (Air FCA)	_	Spusťte zkoušku těsnosti ramena Air FCA a autotest funkce cLLD.
Air FCA MultiSense	Proveďte autotest funkce cLLD ramena Air FCA MultiSense	_	Spusťte zkoušku těsnosti ramena Air FCA a autotest funkce cLLD.



Přístroj / Komponenta	Úloha péče o systém	Čisticí prostředek / Jednorázový výrobek / Zařízení	Reference / Činnosti péče o systém
Nádoba na systémovou kapalinu	Vyčistěte	Deionizovaná voda, alkohol, slabý čisticí prostředek, povrchově aktivní látka, dezinfekční prostředek, zásada, bělidlo	Viz část "Čištění nádoby systémové kapaliny a nádoby na odpad" [▶ 142].
Mycí stanice (Liquid FCA)	Vyčistěte	Detergent nebo antiseptický roztok	_
Podložky prstů upínadla RGA	Odstraňte částice a zbytky z podložek prstů upínadla	Hadřík bez žmolků s alkoholem	Otřete čisticím prostředkem.
Dokovací stanice a prsty upínadla (rozhraní pro připevnění)	Odstraňte částice a zbytky z rozhraní pro upevnění prstů upínadla (PCBA, magnet a kužel)	Hadřík bez žmolků s alkoholem	Otřete čisticím prostředkem.
Okno samostatného skeneru čárových kódů	Vyčistěte	Slabý detergent	VAROVÁNÍ! Nedívejte se do laserového paprsku. Viz návod výrobce skeneru čárových kódů. Viz část "Přístroj s laserovým zářením" [▶ 38].
Okno skeneru Fluent ID a rotátoru zkumavek	Zkontrolujte, zda není znečištěné a poškozené V případě potřeby vyčistěte	Slabý detergent Deionizovaná voda pro oplach	VAROVÁNÍ! Nedívejte se do laserového paprsku. Vyčistěte a opláchněte pomocí měkkého hadříku.
Reflektor Fluent ID a rotátoru zkumavek	Zkontrolujte, zda není znečištěné a poškozené V případě potřeby vyčistěte	Slabý detergent Deionizovaná voda pro oplach	VAROVÁNÍ! Nedívejte se do laserového paprsku. Vyčistěte a opláchněte pomocí měkkého hadříku.
Upínadlo FCA	Vyčistěte	Alkohol	_
Rotátor zkumavek	Zkontrolujte přítomnost a těsnost blokovacích a polohovacích kolíků. V případě potřeby kolíky dotáhněte nebo vyměňte.	_	Viz část "Výměna blokovacích kolíků a polohovacích kolíků" [▶ 194].



Přístroj / Komponenta	Úloha péče o systém	Čisticí prostředek / Jednorázový výrobek / Zařízení	Reference / Činnosti péče o systém
MCA 96	Proveďte zkoušku těsnosti	-	Spusťte metodu kontroly těsnosti ramena MCA 96 .
MCA 96	Proveďte test výkonu pipetování	_	 Spusťte metodu kontroly výkonu pipetování ramena MCA 96, pokud: Životnost těsnění kužele dosáhla 90 % Jsou používány špičky 10 ul nebo 50 ul
			Po dosažení 90 % životnosti těsnění kužele vydá systém Fluent Control varování.

7.4.3 Měsíční péče o systém

V následující tabulce jsou uvedeny úkoly měsíční péče o systém v chronologickém pořadí:

Tab. 25: Tabulka měsíční péče o systém

Přístroj / Komponenta	Úloha péče o systém	Čisticí prostředek / Jednorázový výrobek / Zařízení	Reference / Činnosti péče o systém
Software	Restartujte počítač	_	Vypněte počítač. Počkejte 10 sekund. Znovu zapněte počítač.
Vodítko ramena	Vyčistěte	Bavlněný tampon nebo hadřík, který nepouští vlákna, na šroubováku	Viz část "Čištění vodítka ramena" [⊁ 148].
MCA 96	Čištění MCH 96	Hadřík bez žmolků s alkoholem, stlačený vzduch	Viz část "Čištění MCH 96" [▶ 150].

7.4.4 Pravidelná péče o systém



Intervaly, ve kterých je třeba tyto úkoly provádět, by měl určit hlavní operátor.

V následující tabulce jsou uvedeny úkoly péče o systém v chronologickém pořadí:



Tab. 26: Tabulka pravidelné péče o systém

Přístroj / Komponenta	Úloha péče o systém	Čisticí prostředek / Jednorázový výrobek / Zařízení	Reference / Činnosti péče o systém
Připojení s kuželovou objímkou	Odstranění částic Vyčištění povrchu	Alkohol, hadřík bez žmolků	
Světlo UVC	Zkontrolujte otisky prstů. V případě potřeby vyčistěte.	Alkohol, hadřík bez žmolků	
Prsty upínadla ramena MCA 96	Vizuální kontrola prstů upínadla zejména poté, co byly vystaveny nárazu	Zkontrolujte, zda nejsou prsty upínadla znečištěné nebo poškozené. Pokud jsou znečištěné, vyčistěte je alkoholem a hadříkem, který nepouští vlákna.	

7.4.5 Roční péče o systém

Roční péče o systém pomáhá udržovat přesnost a preciznost a minimalizovat prostoje přístroje. Pomáhá také prodlužovat životnost produktu Fluent.

Obraťte se na místní servisní organizaci Tecan a naplánujte si schůzku pro roční péči o systém. Viz část "Zákaznická podpora" [▶ 199].

7.4.6 Dvouletá péče o systém

Následující úkony preventivní údržby se musí provádět každé 2 roky:

Komponenta	Úkol	Reference
Upínadlo FCA	Vyměňte prsty upínadla FCA. Resetujte čítač ve FluentControl.	Informace o objednávání naleznete v Referenční příručce. Viz "Referenční dokumenty" [> 9].

7.5 Činnosti péče o systém

Chcete-li provést níže popsané činnosti péče o systém, postupujte takto:

- Přepněte do režimu péče o systém. Viz část "Režim péče o systém" [▶ 112].
- Postupujte podle níže uvedených pokynů.



7.5.1 Přemístění přístroje na skříni v rámci laboratoře

▲ UPOZORNĚNÍ

Poškození kabinetu!

Police kabinetu mohly být odstraněny, například kvůli instalaci odstředivky nebo odpadního vozíku. Přemisťování přístroje na kabinetu bez nainstalovaných polic kabinetu může kabinet poškodit a způsobit zranění.

- Před přemístěním přístroje nainstalujte police kabinetu.
- Kabinet přemisťujte pouze po rovné podlaze bez schodů a vyjetých kolejí. Pokud se vyskytují schody nebo vyjeté koleje, použijte zvedací tyče Fluent ke zvednutí systému přes překážku nebo se obraťte na servisního zástupce.

Chcete-li přístroj přemístit na skříni v rámci místnosti, postupujte následovně:

- 1. Ujistěte se, že je skříň bezpečně zaparkovaná a zajištěná proti odjetí.
- 2. Zkontrolujte, zda jsou nainstalovány skříňové police (A).





- 3. Otočte maticí na nohách skříně (B) pomocí klíče.
- 4. Otáčejte červeným šroubem (D) na nohách skříně (C), dokud se zámek neuvolní a kolečka nebudou v pohyblivé poloze.



- 5. Přesuňte přístroj na skříni na nové místo.
- 6. Ujistěte se, že je skříň bezpečně zaparkovaná a zajištěná proti odjetí.



7.5.1.1 Vyrovnání přístroje

Chcete-li přístroj vyrovnat, postupujte následovně:

1. Pomocí klíče spusťte dolů všechny nastavitelné nohy, dokud nebude možné ručně otáčet kolečky skříně.



2. Povolte pojistnou matici (A) na příslušné noze.



 Umístěte referenční segment podle níže uvedených pozic mřížky. Velikost nástroje 480: Pozice mřížky na levé straně 1 a pozice mřížky na pravé straně 21.

Velikost nástroje 780: Pozice mřížky na levé straně 1 a pozice mřížky na pravé straně 41.







Velikost nástroje 1080: Pozice mřížky na levé straně 1 a pozice mřížky na pravé straně 59.

- 4. Pomocí vodováhy se ujistěte, že je přístroj vodorovně a svisle vyrovnán.
- 5. Podle potřeby nastavte úroveň skříně (ve směru hodinových ručiček pro zvýšení, proti směru hodinových ručiček pro snížení).



- 6. Po vyrovnání přístroje utáhněte pojistné matice na nohách skříně.
- 7. Ujistěte se, že je skříň bezpečně zaparkovaná a zajištěná proti odjetí.

7.5.2 Čištění jednorázových kuželů hrotů

Při čištění jednorázového kuželu hrotu postupujte následovně:

- 1. Jednorázové kužely hrotů očistěte alkoholem pomocí hadříku, který nepouští vlákna.
- 2. Během péče o systém zkontrolujte kužely jednorázových hrotů a vyčnívající hrot.

Pro FCA s kapalinou: Ujistěte se, že nástavec prodloužení trubky vyčnívající mimo kužel není poškozen.

3. Zkontrolujte, zda jsou nástavce trubek čisté a bez usazenin.



7.5.3 Čištění fixních hrotů

A UPOZORNĚNÍ

Nebezpečí poranění o fixní hroty při čištění

Pipetovací fixní hroty mohou způsobit poranění.

• Při přístupu k pracovnímu stolu se vyhněte kontaktu s pipetovacími hroty a kontaktu s aerosoly tím, že budete nosit vhodný ochranný oděv.

Při čištění fixních hrotů postupujte následovně:

- 1. Fixní hroty očistěte alkoholem pomocí hadříku, který nepouští vlákna.
- 2. Zkontrolujte, zda jsou fixní hroty čisté a bez usazenin.

7.5.4 Čištění propichovacích hrotů

Chcete-li vyčistit propichovací hroty, spusťte metodu **Údržba čištění propichovacích hrotů**. Tuto metodu je třeba upravit podle nastavení vašeho pracovního stolu.

Skript obsahuje následující kroky:

- 1. Připravte pracovní stůl (tj. laboratorní vybavení a hardware).
- 2. Propíchněte do Z-startu 8 prázdných zkumavek s uzávěrem na rotátoru zkumavek nebo na držáku zkumavek.
- Přístupnou část propichovacích hrotů ručně očistěte 70% etanolem nebo 2% bělidlem pomocí hadříku, který nepouští vlákna. Vyvarujte se kontaktu s ostrým vrcholem propichovacích hrotů.
- 4. Po ručním čištění proveďte mycí příkazy.

7.5.5 Čištění rotátoru zkumavek

Obecný postup čištění

- K čištění částí rotátoru zkumavek použijte hadříky, které nepouštějí vlákna, a namočte je do jedné z následujících čisticích kapalin: 2% bělidlo, 70% ethanol, 100% isopropanol
- Otřete díly namočenými utěrkami za účelem vyčištění a dezinfekce. K čištění míst, na která nelze dosáhnout hadříkem bez chloupků, použijte vatové tampony.
- 3. Do 5 minut po použití čisticích kapalin je otřete hadříkem namočeným ve vodě.



Demontáž a čištění desky držáku zkumavky

1. Chcete-li uvolnit desku držáku (A), přidržte ji jednou rukou a druhou rukou zatáhněte za blokovací kolík držáku (B).

2. Vyjměte desku držáku z rotátoru zkumavek.



 Vyčistěte desku držáku podle výše uvedených obecných pokynů nebo můžete držák zkumavky inkubovat v lázni s 2% bělidlem po dobu maximálně 2 hodin.

Čištění povrchů rotátoru zkumavek

1. Vyčistěte přístupné povrchy rotátoru zkumavek podle výše uvedených obecných pokynů.





2. Chcete-li změnit polohu bubnu ručně, přidržte buben jednou rukou a stiskněte uvolňovací tlačítko solenoidu.

- 3. Otáčejte bubnem ručně a uvolněte uvolňovací tlačítko solenoidu.
- 4. Otáčejte bubnem, dokud jej solenoid nezablokuje.
- 5. Vyčistěte povrchy, které předtím nebyly přístupné, podle výše uvedených obecných pokynů.

Montáž desky držáku

- 1. Umístěte desku držáku na horní část bubnu rotátoru zkumavek.
- 2. Jednou rukou přitlačte desku držáku ke spodní části přístroje a zatlačte černý jezdec dozadu, aby se deska držáku zajistila.





7.5.6 Čištění mycí stanice rotátoru zkumavek

Obecný postup čištění

- Mycí stanici lze čistit na pracovním stole nebo ji lze pro účely čištění demontovat.
- ✓ Pro lepší přístup používejte místo hadříků kartáč na láhve.
- K čištění částí mycí stanice použijte hadříky, které nepouštějí vlákna, a namočte je do jedné z následujících čisticích kapalin: 2% bělidlo, 70% ethanol, 100% isopropanol
- 2. Otřete díly namočenými utěrkami za účelem vyčištění a dezinfekce.
- 3. Do 5 minut po použití čisticích kapalin je otřete hadříkem namočeným ve vodě.

Demontáž mycí stanice rotátoru zkumavek

1. Stiskněte uvolňovací páčku mycí stanice směrem ke skříni skeneru čárových kódů a druhou rukou mycí stanici zvedněte.







2. Odpojte odpadní trubky a vložte konektory do držáků odpadních trubek.

Montáž mycí stanice rotátoru zkumavek

1. Připojte konektory odpadních trubek.









 Namontujte mycí stanici na vodicí hřídele (A) a přitlačte ji na základní desku. Zkontrolujte, zda uvolňovací páčka zapadla zpět na místo a drží mycí stanici v poloze.

7.5.7 Čištění podložek a segmentů

Při čištění podložek a segmentů postupujte následovně:

OZNÁMENÍ

Selhání detekce kapaliny (cLLD)!

Možná porucha detekce kapaliny (cLLD) v důsledku narušeného kontaktu mezi podložkou a segmentem plošiny.

Vždy se ujistěte, že jsou podložky a segmenty čisté a suché.

- Vyjměte podložky z plošiny přístroje. Segmenty a soupravy se čistí na místě.
- 2. Otřete povrch podložek, segmentů a souprav čisticím prostředkem.

Podložky, segmenty a soupravy opláchněte DI vodou.

3. Vraťte podložky zpět na plošinu přístroje.

7.5.8 Čištění podnosů plošiny

Při čištění podnosů plošiny postupujte následovně:

- ✓ Segmenty nad podnosem plošiny jsou demontovány. Viz "Odebírání segmentů" [▶ 83].
- Pokud segmenty plošiny, jako je Fluent ID nebo aktivní nosič MCA, nelze demontovat, posuňte podnosy plošiny do otevřené polohy.
- 1. Demontujte podnosy plošiny z přístroje.



- 2. Vyprázdněte podnosy odstraněním kapaliny v souladu s laboratorním protokolem pro manipulaci s danou kapalinou.
- 3. Pokud dojde k poškození nebo ztrátě podnosů plošiny, je třeba je vyměnit.
- 4. Otřete povrch podnosů plošiny čisticím prostředkem.
- Vložte podnosy plošiny zpět do přístroje.
 Natočte podnosy plošiny podle obrázku níže.
 Sousední podnosy plošiny musí být vzájemně propojeny.



Obr. 44: Nesprávné umístění podnosu plošiny



Obr. 45: Zablokování podnosů plošiny





Obr. 46: Správné umístění podnosu plošiny

7.5.9 Výměna reflexní fólie Fluent ID

- ✓ Samolepicí reflexní fólie
- 1. Zahřejte reflexní fólii. Použijte horkovzdušnou pistoli.
- 2. Sejměte reflexní fólii.



3. Případné zbytky odstraňte alkoholem.





4. Na horní konec reflektoru nalepte novou samolepicí reflexní fólii.

7.5.10 Aplikace reflexní fólie na odpadní skluz DiTi

- ✓ Samolepicí reflexní fólie
- 1. Přiložte pásku na odpadní skluz DiTi podle obrázku níže.



2. Na odpadní skluz DiTi nalepte novou samolepicí reflexní fólii podle obrázku níže.





Laserový paprsek musí být uprostřed reflexní fólie.

3. Odstraňte pásku z odpadního skluzu DiTi.

7.5.11 Čištění bezpečnostních panelů

Při čištění bezpečnostních panelů postupujte následovně:

1. Otřete vnitřní a vnější povrch bezpečnostních panelů čisticím prostředkem.

7.5.12 Čištění jednotky odpadu z jednorázových špiček a mycí stanice

Chcete-li vyčistit odpad z jednorázových hrotů a jednotku mycí stanice, postupujte následovně:

- 1. Stiskněte tlačítko rychlovypínacího uzávěru (B).
- 2. Posuňte mycí stanici dozadu.



Obr. 47: Upevňovací prvek pro vyjmutí pouzdra sáčku

3. Demontujte mycí stanici z odpadu jednorázových špiček a z jednotky mycí stanice.





Obr. 48: Demontáž mycí stanice

- 4. Otřete povrch mycí stanice čisticím prostředkem a odstraňte rozlitá činidla.
- 5. Stiskněte tlačítko rychlovypínacího uzávěru (B).
- 6. Umístěte mycí stanici (A) na místo.
- 7. Zatlačte mycí stanici dopředu.



Obr. 49: Zpětná montáž mycí stanice

7.5.13 Čištění skluzu odpadních jednorázových hrotů

Při čištění skluzu odpadních jednorázových hrotů postupujte následovně:

- ✓ Přední bezpečnostní panel je otevřený.
- 1. Sejměte kryt (A) ze skluzu odpadních jednorázových hrotů.





Obr. 50: Demontáž krytu ze skluzu odpadních jednorázových hrotů

2. Sejměte skluz odpadních jednorázových hrotů (B) z držáku.



Obr. 51: Demontáž skluzu odpadních jednorázových hrotů

3. Přidržte utěrku pod spodním otvorem (C) skluzu odpadních jednorázových hrotů.

Zabraňte odkapávání kontaminovaných látek.



Obr. 52: Manipulace se skluzem odpadních jednorázových hrotů

- 4. Umístěte skluz odpadních jednorázových hrotů a kryt do nádoby naplněné čisticím prostředkem.
- 5. Nechte je namočené 30 minut až 4 hodiny.
- 6. Vyjměte z nádoby skluz odpadních jednorázových hrotů a kryt a položte je na čistý suchý ručník.



- 7. Nechte uschnout.
- 8. Namontujte zpět skluz odpadních jednorázových hrotů (B) na držák.



- Obr. 53: Nainstalujte skluz odpadních jednorázových hrotů zpět do sady
- 9. Zkontrolujte, zda je polohovací kolík správně zasunutý do drážky (D).
- 10. Umístěte kryt (A) na horní část odpadního skluzu.



Obr. 54: Polohovací kolík a kryt

7.5.14 Výměna sáčku na odpad jednorázových hrotů

Při výměně sáčku odpadních jednorázových špiček postupujte následovně:

1. Zvedněte upevňovací prvek (A) a posuňte kryt sáčku dopředu.



Obr. 55: Upevňovací prvek pro vyjmutí pouzdra sáčku



- 2. Vyjměte kryt jednorázového sáčku (A).
- 3. Vyjměte jednorázový odpadní sáček na špičky (B).
- 4. Zlikvidujte jednorázový odpadní sáček na špičky podle pokynů vaší laboratoře.
- 5. Vložte nový jednorázový sáček na odpad (B) do prázdného pouzdra sáčku (B).



Obr. 56: Pouzdro na sáčky a jednorázový sáček na odpad



Obr. 57: Správná montáž odpadního skluzu na pracovním stole



A UPOZORNĚNÍ

Nesprávně umístěný odpadní skluz může způsobit náraz ramena nebo nesprávné vysunutí špiček. Odpadní skluzy musí být správně umístěny, jak je znázorněno níže: vložte obrázky správného/nesprávného umístění odpadních skluzů

6. Zasuňte kryt sáčku na místo a uzavřete jej pomocí uzávěru (A).





Obr. 58: Uzavírací upínací prvek

7.5.15 Čištění cesty kapaliny

- Chcete-li odstranit zbytky bílkovin z vnitřku fixních hrotů, použijte slabou kyselinu a následně zásaditý čisticí prostředek.
- K odstranění zbytků nukleových kyselin z vnitřku fixních hrotů použijte zásaditý čisticí prostředek.
- Proces mohou ovlivnit čisticí prostředky, například Decon/Contrad. Pokud tedy takové prostředky používáte, pečlivě proces validujte.
- Isopropanol je vysoce účinný dezinfekční prostředek. Rychle se odpařuje a zanechává povrchy připravené k použití.
- Používejte pouze povolené čisticí prostředky. K proplachování celého kapalinového systému nepoužívejte bělicí roztoky.

Při čištění cesty kapaliny postupujte následovně:

- 1. Odpojte trubku systémové kapaliny od nádoby na kapalinu.
- 2. Připojte trubku pro údržbu (30043739) k systémovému potrubí.
- 3. Umístěte otevřený konec trubky pro údržbu do láhve s čisticím prostředkem.
- 4. Propláchněte čisticím prostředkem (20 ml s RapidWash a 10 ml s ředidlem).
- 5. Nechte je namočené 20 minut.
- 6. Vložte trubku do láhve s DI vodou.
- 7. Dvakrát opláchněte DI vodou (20 ml s RapidWash a 10 ml s ředidlem).

A VAROVÁNÍ

Hořlavé kapaliny!

Nebezpečí požáru způsobené hořlavými kapalinami nebo kapalinou v systému.

- Zabraňte vzniku a hromadění hořlavých výparů.
- Neprovozujte systém bez podnosů plošiny.
- 8. Vložte trubku do láhve s alkoholem.
- 9. Propláchněte alkoholem (20 ml s RapidWash a 10 ml s ředidlem).
- 10. Odstraňte trubku pro údržbu ze systémové trubky a připojte systémovou trubku k nádobě systémové kapaliny.



- 11. Dvakrát propláchněte DI vodou (20 ml s přípravkem RapidWash a pětinásobek objemu ředidla).
- 12. Zkontrolujte, zda se v trubce netvoří bubliny.
- 13. Pokud se objeví bublinky, znovu je propláchněte.

7.5.16 Připojení nádoby systémové kapaliny a nádoby na odpad

Při přípravě nádoby systémové kapaliny a nádoby na odpad postupujte následovně:

6

Bezproblémový provoz je zaručen pouze v případě použití originálních nádob s řídicím systémem od společnosti Tecan. Před prvním použitím je třeba nádobu se systémovou kapalinou důkladně ručně vypláchnout, aby se z vnitřku láhve odstranily veškeré pevné zbytky. Viz část "Čištění nádoby systémové kapaliny a nádoby na odpad" [▶ 142].

- ✓ Nádoba Tecan o objemu více než 20 litrů
- 1. Zkontrolujte, zda je systém detekce kapalin (A, B) správně připojen.
- 2. Zkontrolujte, zda jsou trubky (C, D) správně připojeny.



7.5.17 Čištění nádoby systémové kapaliny a nádoby na odpad

Při čištění nádoby kapaliny a nádoby na odpad postupujte následovně:

- 1. Vyprázdněte ručně nádobu na mycí kapalinu.
- 2. Vyčistěte nádobu na kapalinu v umyvadle čisticím prostředkem a opláchněte ji.
- 3. Nádobu na kapalinu vydezinfikujte alkoholem.
- Připojte nádobu na systémovou kapalinu a nádobu na odpad, viz část "Připojení nádoby systémové kapaliny a nádoby na odpad" [▶ 142].

7.5.18 Kontrola těsnosti stříkaček

Při kontrole správné těsnosti stříkaček postupujte následovně:





Obr. 59: Kontrola těsnosti

A Pojistný šroub pístu

B Šroub stříkačky

- Posuňte písty do středu stříkaček pomocí skriptu údržby, který nasává vzduch. Poznámka: Skript údržby vám poskytne správce systému FluentControl laboratoře.
- 2. Utáhněte šroub stříkačky (B), tj. otočte jej doprava.
- 3. Otáčením pojistného šroubu pístu (A) ve směru hodinových ručiček jej utáhněte.

7.5.19 Kontrola těsnění (MCA 384)

Při kontrole těsnění (MCA) postupujte takto:

1. Zkontrolujte, zda těsnění (A) není poškozené.



Viz část "Výměna těsnění (MCA 384)" [> 144].

Ujistěte se, že těsnění nejsou poškozená. Poškozené těsnění je třeba vyměnit.



7.5.20 Výměna těsnění (MCA 384)



В

D

Obr. 60: Díly a nástroje

- A Těsnění
- **C** Nástroj na vložení tupé trubky
- E Nástroj na vyjmutí tupé trubky

Při výměně těsnění (MCA) postupujte takto:

- 1. Sejměte desku adaptéru.
- 2. Přesuňte pipetovací hlavu dopředu a zvedněte ji co nejvýše.
- 3. Zatlačte nástroj na vyjmutí tupé trubky asi 2 mm do kanálu.
- Tupá trubka

Nástroj na vložení těsnění




4. Tupou trubku zajistěte otáčením nástroje ve směru hodinových ručiček.

5. Vytáhněte tupou trubku z kanálu.









6. Pomocí dlouhých kleští vyjměte těsnění z kanálu.

- 7. Nové těsnění namažte vrstvou minerálního oleje.
- 8. Pomocí nástroje pro vkládání těsnění vložte namazané těsnění do kanálu.



9. Vnější povrch nové tupé trubky namažte minerálním olejem.





10. Pomocí nástroje pro vkládání tupých trubek opatrně zasuňte namazanou tupou trubku do kanálu.

Tupá trubka musí být zcela zasunuta do kanálu.



- 11. Zkontrolujte, zda je nové těsnění vyrovnané s ostatními těsněními.
- 12. Spusťte metodu Zkouška těsnosti MCA.

7.5.21 Čištění bloku fixních špiček (MCA 384)

- Blok fixních špiček (MCA) očistěte alkoholem pomocí hadříku, který nepouští vlákna.
- 2. Blok špiček otřete do sucha utěrkou, která nepouští vlákna, nebo blok špiček vysušte stlačeným vzduchem bez obsahu oleje.
- 3. Zkontrolujte, zda je blok fixních špiček (MCA) čistý a bez usazenin.
- 4. Blok špiček uložte do krabice na bloky špiček.

Aby se zabránilo kontaminaci při manipulaci s blokem špiček:

- Blok špiček musí být uložen na bezprašném místě.
- Nikdy se špiček nedotýkejte prsty. Blok špiček musí být při manipulaci vždy přidržován blokem PEEK.
- Blok špiček nikdy nepokládejte špičkami na stůl.

7.5.22 Čištění adaptéru desky (MCA 384)

Při čištění adaptéru desky postupujte následovně:

- 1. Zkontrolujte, zda nejsou v kanálech nečistoty.
- 2. K odstranění případné mastnoty použijte alkohol.
- 3. Vysušte stlačeným vzduchem.



7.5.23 Čištění vodítka ramena

Při čištění vodítka ramena postupujte následovně:

- 1. Vyčistěte váleček vodítka ramena (A) na vodítka ramena vatovým tamponem nebo hadříkem bez žmolků na šroubováku.
- 2. Vyčistěte kolejnice ramena (B) hadříkem, který nepouští vlákna.
- 3. Horní povrch vodicí lišty na vodítku ramena MCA očistěte hadříkem, který nepouští vlákna.



7.5.24 Utahování kužele DiTi

Při utahování kužele DiTi ramena FCA postupujte následovně:

- 1. Přidržte adaptér hrotu (D) a trubku pro vyhazování hrotu (C).
- 2. Kužel DiTi (A) utáhněte pomocí klíče na kužel DiTi (B).



3. Spusťte metodu Rutinní údržba FCA .



7.5.25 Čtečka Frida

Vložka



Obr. 61: Vložka čtečky Frida Při instalaci nasaďte vložku do čtečky Frida a vyrovnejte značky.

Slepá zátka



Obr. 62: Slepá zátka čtečky Frida

Slepá zátka chrání čtečku Frida při vyjmutí vložky. Pro instalaci zasuňte slepou zátku do čtečky Frida.



7.5.26 Čištění MCH 96

Zkontrolujte ejekční desku (A) ohledně viditelného poškození. Rozlití určitých provozních kapalin (například DMSO nebo acetonitrilu) nebo čisticích prostředků (například bělicích prostředků) může vést k poškození desky. V takovém případě požádejte místní servisní organizaci o výměnu.



- 1. Pomocí nástroje pro přemístění (Move Tool) posuňte ejekční osu do spodní části. 2. . 3.
- K čištění horního povrchu ejekční desky (C), kuželové desky a kuželů (B) použijte stlačený vzduch bez příměsi oleje
- 3. Hadříkem bez žmolků s alkoholem vyčistěte spodní povrch ejekční desky (D).









Obrázek níže vlevo zobrazuje ejekční desku před čištěním a vpravo po čištění.



8 Řešení problémů

V této kapitole naleznete nápovědu k obnovení provozu po problému s přístrojem Fluent. Další informace nebo informace pro případ problémů, které nejsou popsány v tomto návodu nebo nejsou dostatečně podrobné, naleznete v části "Zákaznická podpora" [> 199].

8.1 Bezpečnostní pokyny pro tuto kapitolu

A UPOZORNĚNÍ

Křížová kontaminace v důsledku poškozených hrotů po havárii!

Ohnuté špičky nebo poškozený povlak špiček způsobují nepřesnost pipetování a chyby při detekci kapalin.

• Zkontrolujte fixní hroty po havárii. Viz část "Kontrola fixních hrotů" [▶ 175].

8.2 Tabulky pro řešení problémů

V tabulkách pro řešení problémů jsou uvedeny možné problémy, jejich příčiny a nápravná opatření. Další informace nebo informace pro případ problémů, které nejsou popsány v této příručce nebo nejsou dostatečně podrobné, naleznete v části "Zákaznická podpora" [> 199].

8.2.1 Řešení problémů s přístrojem

Tab. 28: Tabulka pro řešení problémů s přístrojem

Problém/chyba	Možná příčina	Nápravné opatření
Úniky systémové kapaliny	Trubky a/nebo spoje trubek jsou netěsné. Injekční stříkačka je netěsná.	Viz část "Zákaznická podpora" [▶ 199].
Chyba komunikace	Napájení není zapnuté. Je přerušeno napájení nebo komunikace. Žádná komunikace.	Vypněte přístroj. Počkejte, až zhasne kontrolka stavu přístroje a kontrolka napájení. Vypněte počítač. Zkontrolujte kabel a zástrčky. Zapněte přístroj a počítač.
	Zablokování pohonu X, Y nebo Z.	Zkontrolujte, zda nejsou překážky. OZNÁMENÍ Ujistěte se, že se ramena mohou volně pohybovat.



Problém/chyba	Možná příčina	Nápravné opatření
Chyba inicializace	Ramena nelze inicializovat.	Zkontrolujte, zda nejsou překážky. OZNÁMENÍ Ujistěte se, že se ramena mohou volně pohybovat.
	Problém hardwaru.	Viz část "Zákaznická podpora" [▶ 199].
Snímač předního bezpečnostního panelu a zámek dveří jsou poškozené	Mechanická porucha zámků dveří.	Vypněte přístroj. Viz část "Zákaznická podpora" [▶ 199].
Chybějící nebo poškozený bezpečnostní panel	Bezpečnost nelze zaručit.	Vypněte přístroj. Viz část "Zákaznická podpora" [▶ 199].
Porucha detekce kapaliny (cLLD)	Znečištěný kontaktní povrch. Nedostatečný kontakt mezi laboratorním vybavením a segmentem.	Připravte plošinu. Viz část . Vyčistěte kontaktní plochu. Viz část "Čištění podložek a segmentů" [▶ 132].
	Systémová kapalina má vodivost > 10 µS/ cm pro kompatibilitu s cLLD.	Obraťte se na hlavního operátora.

8.2.2 Řešení problémů s flexibilním ramenem kanálu (FCA)

Tab. 29: Tabulka řešení problémů s flexibilním pipetovacím ramenem

Problém/chyba	Možná příčina	Nápravné opatření
Uvolněný kužel DiTi UPOZORNĚNÍ! Nepřesné pipetování objemů!	Nedostatečně utažený kužel DiTi.	Utáhněte kužel DiTi.
Jednorázová špička není vyzvednuta	Nedostatečně utažený kužel DiTi.	Utáhněte kužel DiTi.



Problém/chyba	Možná příčina	Nápravné opatření
Jednorázová špička není zlikvidována	Nedostatečně utažený kužel DiTi.	Utáhněte kužel DiTi.
	Znovu použité DiTi	Ujistěte se, že jsou špičky nové. DiTi se nedoporučuje používat opakovaně.
		Jednorázové špičky nevyhozeny do odpadního skluzu / odpadní skluz není správně umístěn
Jednorázové špičky nevyhozeny do odpadního skluzu	Odpadní skluz není správně umístěn	Zajistěte správné umístění odpadních skluzů. Viz část "Čištění skluzu odpadních jednorázových hrotů" [▶ 137]
Špičky nejsou zarovnány	Nosič ve špatné poloze.	Zajistěte správnou polohu nosiče. Viz část "Vkládání standardních podložek" [▶ 85].
s laboratorním vybavením na jednom nosiči	Segment není uzamčen na místě. Laboratorní vybavení není správně umístěno.	Uzamkněte segment na místě. Viz část "Kontrola segmentu" [▶ 192].
Špičky nejsou zarovnány s laboratorním vybavením na několika nosičích	Vadné seřízení ramena způsobené kolizí.	Viz část "Zákaznická podpora" [▶ 199].
Špička naráží na dno laboratorního vybavení	Nesprávné laboratorní vybavení. Laboratorní vybavení není správně umístěno.	Ujistěte se, že laboratorní vybavení na pracovní ploše odpovídá uspořádání pracovní plochy dané metody.
DiTi odkapává	Zašpiněný kužel DiTi způsobuje netěsnost.	Vyčistěte kužel DiTi.
	Znovu použité DiTi	Ujistěte se, že jsou špičky nové. DiTi se nedoporučuje používat opakovaně.
Chybové hlášení: Tlak mimo rozsah (Air	Mokrý inline filtr po aspiraci s nesprávnou velikostí DiTi.	Ujistěte se, že velikost DiTi na plošině odpovídá velikosti definované v metodě. Ujistěte se, že jsou kužely DiTi správně utaženy
FCA)		Zkontrolujte inline filtr. Viz část "Kontrola řadového filtru (FCA se vzduchem)" [▶ 168].



Problém/chyba	Možná příčina	Nápravné opatření
Chybová hlášení:	Magnetické pole ruší snímač přítomnosti	_
DiTi nepřineseno	DiTi.	
DiTi neupuštěno		
Detekce kapaliny: Špička nedetekuje kapalinu	Uvolněný kužel DiTi	Utáhněte kužel DiTi (viz část "Utahování kužele DiTi" [▶ 148]).
Opotřebované O-kroužky	Adaptéry špiček MultiSense	Vyměňte O-kroužky a X-kroužek na adaptérech špiček MultiSense.

8.2.3 Míchání a propichování

Tab. 30: Řešení problémů

Příznak	Možná příčina	Nápravná opatření
Propichovací hrot nelze zatáhnout pomocí softwarových příkazů	Zaseknutý propichovací hrot	Viz část "Vytahování zaseknutých propichovacích hrotů" [▶ 186].
Poškozený propichovací hrot	Ohnutý propichovací hrot Poškozený hrot	Vyměňte propichovací hrot. Viz části "Demontáž propichovacích hrotů" [▶ 179] a "Montáž propichovacích hrotů" [▶ 182].



Příznak	Možná příčina	Nápravná opatření
Chyby propichování	Propichovací hrot je příliš suchý	Navlhčení vodou (mycí stanice)
	Poškozený hrot	Vyměňte propichovací hrot.
	Ohnutý propichovací hrot	VIZ casti "Demontaz propichovacích hrotů" [▶ 179] a "Montáž propichovacích hrotů" [▶ 182].
	Špatný parametr propichování	Obraťte se na hlavního operátora.
	Použití nesprávného typu pohybu	Obraťte se na hlavního operátora.
	Použití nesprávných zkumavek	Používejte podporované zkumavky. Viz část "Podložky rotátoru zkumavek" [▶ 63].
	Rameno dosáhlo své životnosti	Viz část "Zákaznická podpora" [▶ 199].
Problémy při manipulaci s kapalinami	Ucpané propichovací hroty	Propláchněte propichovací hroty. Obecně zkontrolujte postup mytí.
	Poškozený hrot	Vyměňte propichovací hrot. Viz části "Demontáž propichovacích hrotů" [> 179] a "Montáž propichovacích hrotů" [> 182].
	Nesprávně namontované injekční stříkačky	Zkontrolujte těsnost stříkaček. Viz část "Kontrola těsnosti stříkaček" [▶ 142].
	Bubliny v kapalinovém systému	Propláchněte. Viz část "Zákaznická podpora" [▶ 199].



Příznak	Možná příčina	Nápravná opatření
Problémy s hemolýzou	Ředění vzorku	Vyšší přebytečný objem nebo rozdělovací objem
		Fyziologický roztok 0,9 % jako rozdělovací objem
		Nižší rychlost pipetování
	Poškozený hrot	Vyměňte propichovací hrot. Viz části "Demontáž propichovacích hrotů" [▶ 179] a "Montáž propichovacích hrotů" [▶ 182].
	Parametry míchání	Ujistěte se, že použité parametry rotace/oscilace rotátoru zkumavek nevedou k hemolýze



Příznak	Možná příčina	Nápravná opatření
Vzorek ve stříkačce	Jakýkoliv	Vyčistěte systém. Viz část "Čištění cesty kapaliny" [▶ 141].
	Špatná vzduchová mezera.	Ověřte postup mytí.
	Nesprávně namontované injekční stříkačky.	Zkontrolujte těsnost stříkaček. Viz část "Kontrola těsnosti stříkaček" [▶ 142]
		Spusťte metodu Metoda úniku FCA.
		Větší přední vzduchová mezera.
		Nižší rychlost aspirace.
	Nesprávná příprava vzorku pro manipulaci s kapalinou. Zkumavky se zdroji vzorků obsahují pevné částice, jako jsou sraženiny, zbytky buněk atd.	Zajistěte správnou přípravu vzorku, aby bylo možné pipetovat kapalinu vzorku.
		Ujistěte se, že zkumavky se zdrojem vzorku neobsahují pevné částice, jako jsou sraženiny, zbytky buněk atd.
	Nesprávná příprava vzorku pro manipulaci s kapalinou. Zkumavky nejsou správně naplněny a stále obsahují částečný podtlak, který snižuje vedoucí vzduchovou mezeru při propíchnutí.	Ujistěte se, že jsou zkumavky se zdroji vzorků správně naplněny cílovým objemem zkumavky.
		Ujistěte se, že zkumavky se zdroji vzorků neobsahují vakuum.
		Zvětšete přední vzduchovou mezeru, abyste vyrovnali případný zbývající podtlak.
Propichovací hrot se v postupu mytí ohýbá	Propichovací hrot není vycentrován v otvorech čističe mycí stanice	Vytvořte kopii promývací stanice a naučte pozice pipetování.



Příznak	Možná příčina	Nápravná opatření
Nesprávná detekce hladiny kapaliny: pouze v určitých kanálech	Ohnutý propichovací hrot: Propichovací hrot je ohnutý, a proto se při propichování dotýká stěny zkumavky	Vyměňte propichovací hrot. Viz části "Demontáž propichovacích hrotů" [▶ 179] a "Montáž propichovacích hrotů" [▶ 182].
	Poloha propichování je nesprávná, a proto se propichovací hrot při propichování dotýká stěny zkumavky.	Používejte laboratorní vybavení vyrobené společností Tecan. Viz část "Podložky rotátoru zkumavek" [▶ 63].
		Výuka/úprava polohy pipetování laboratorního vybavení
	Orientace propichovacího hrotu je nesprávná.	Nasaďte propichovací hrot otvorem směrem k přední části přístroje. Viz část "Montáž propichovacích hrotů" [▶ 182].
Nesprávná detekce hladiny kapaliny: stálá odchylka očekávané a detekované hladiny kapaliny	Výrobní tolerance ramena, rotátoru zkumavek a/nebo přístroje v kombinaci s propichovací silou mohou vést ke znatelnému posunu Z při detekci hladiny kapaliny.	Naučte/upravte vlastní atribut "PiercingDetectionHeightCom pensation" v definici laboratorního vybavení zkumavky

8.2.4 Řešení problémů s vícekanálovým ramenem (MCA)

Tab. 31: Tabulka řešení problémů s vícekanálovým ramenem kanálu

Problém/chyba	Možná příčina	Nápravné opatření
Hroty nejsou	Mechanická závada	Viz část "Zákaznická podpora" [▶ 199].
vyrovnany s nosici	Havárie ramena	
Mikrodeska a pipetovací hlava nejsou 100% paralelní	Srážka	Obraťte se na hlavního operátora, aby zkontroloval rovnoběžnost pipetovací hlavy a plošiny.
Během pipetování přestane pipetovací hlava generovat chybu	Zrychlení aspirace a dávkování je v porovnání s rychlostí příliš rychlé.	Zrychlení musí být v přiměřeném poměru k rychlosti aspirace a dávkování.
	Zpomalení aspirace a dávkování je v porovnání s rychlostí příliš rychlé.	Zpomalení musí být v přiměřeném poměru k rychlosti aspirace a dávkování.
		Problém nelze vyřešit. Viz část "Zákaznická podpora" [▶ 199].



Problém/chyba	Možná příčina	Nápravné opatření
Netěsnost několika nebo všech pipetovacích kanálů	Špatné jednorázové hroty, blok fixních hrotů, adaptér fixního hrotu, těsnění kuželu hrotu nebo těsnění.	Vždy používejte jednorázové hroty, blok fixních hrotů, adaptér fixních hrotů, těsnění kuželů hrotů nebo těsnění dodávané společností Tecan. Obraťte se na hlavního operátora, aby systém propláchl a zkontroloval jeho těsnost.
	Těsnění kuželu hrotu nebo další těsnění jsou stará nebo vadná.	Viz část "Zákaznická podpora" [▶ 199]. Obraťte se na hlavního operátora, aby systém propláchl a zkontroloval jeho těsnost.
	Pipetovací hlava je vadná.	Viz část "Zákaznická podpora" [▶ 199].
Únik z jednoho kanálu	Těsnění kuželu hrotu nebo jiná těsnění v pipetovací hlavě jsou vadná.	Obraťte se na hlavního operátora, aby systém propláchl a zkontroloval jeho těsnost. Viz část "Zákaznická podpora" [▶ 199].
Jeden jednorázový hrot není správně vyzvednut	Jednotlivý jednorázový hrot je vadný. Těsnění kuželu hrotu na této pozici jednorázového hrotu je vadné.	Vyměňte jednorázové hroty. Problém nelze vyřešit. Viz část "Zákaznická podpora" [▶ 199].
Jeden jednorázový hrot se neupouští	Jednotlivý jednorázový hrot je vadný. Těsnění kuželu hrotu na této pozici jednorázového hrotu je vadné.	Problém nelze vyřešit. Viz část "Zákaznická podpora" [▶ 199].
	Špatná vlhkost	Zkontrolujte, zda je vlhkost v mezích provozní vlhkosti. Viz část "Podmínky prostředí" [> 44].
Několik nebo všechny jednorázové hroty se neupouští	Použití nesprávných jednorázových hrotů.	Vždy používejte jednorázové hroty dodávané společností Tecan. Problém nelze vyřešit. Viz část "Zákaznická podpora" [▶ 199].



Problém/chyba	Možná příčina	Nápravné opatření
Při vyzvedávání jednorázových hrotů se box s jednorázovými hroty zvedne	Nosič není správně nastaven. Chybně zadaný posun X a/nebo Y.	Nastavte přesně všechny nosiče (mechanické). Vyměňte nosič jednorázových hrotů. Problém nelze vyřešit. Viz část "Zákaznická podpora" [▶ 199].
	Box jednorázových hrotů nesplňuje specifikace.	Vždy používejte boxy jednorázových hrotů, které odpovídají standardům Společnosti pro biomolekulární screening. Problém nelze vyřešit. Viz část "Zákaznická podpora" [▶ 199].
	Nosič jednorázových hrotů je vadný (závada na držácích jednorázových hrotů).	Vždy používejte boxy jednorázových hrotů, které odpovídají standardům Společnosti pro biomolekulární screening. Problém nelze vyřešit. Viz část "Zákaznická podpora" [▶ 199].
Nepřesné výsledky pipetování	Jednorázové hroty nejsou správně vyzvednuty. Parametry manipulace s kapalinou jsou nesprávné. Nosiče nejsou správně nastaveny. Pipetovací hlava je vadná.	Obraťte se na hlavního operátora, aby zkontroloval skript aplikace a nosiče. Obraťte se na hlavního operátora, aby zkontroloval parametry prostředí a výšku dávkování. Zkontrolujte povrchovou úpravu hrotu. V případě potřeby vyměňte blok hrotů.
	Povlak standardních hrotů je poškozený.	Obraťte se na hlavního operátora, aby zkontroloval skript aplikace a nosiče. Zkontrolujte povrchovou úpravu hrotu. V případě potřeby vyměňte blok hrotů.
Přenesení	Povlak standardních hrotů je poškozený.	Zkontrolujte povlak. Vyměňte blok fixních hrotů nebo adaptér fixního hrotu.
	Nevhodný skript aplikace.	Kontaktujte hlavního operátora, aby upravil nastavení mytí nebo manipulace s kapalinou a použil jiný mycí roztok.
	Mycí kanály jsou ucpané.	Vyčistěte mycí stanici. Problém nelze vyřešit. Viz část "Zákaznická podpora" [▶ 199].





8.2.5 Řešení problémů s robotickým upínacím ramenem (RGA – Robotic Gripper Arm)

Tab. 32: Tabulka řešení problémů s robotickým upínacím ramenem

Problém/chyba	Možná příčina	Nápravné opatření
Mikrodeska nebyla vyzvednuta	Žádné mikrodesky na nosiči. Prsty upínadla nemohou mikrodesku zvednout.	Položte mikrodesku na nosič. Nastavte polohu upínadla. Vyčistěte prsty upínadla RGA.
Neobvyklý hluk při pohybu ramena	Díly jsou poškozené nebo opotřebované.	Viz část "Zákaznická podpora" [▶ 199].
Nesprávně vyrovnané excentrické upínací prsty	Srážka náhradních prstů. Šrouby prstů nejsou dostatečně utažené.	Vyrovnejte prsty excentrického upínadla. Viz část "Kontrola vyrovnání prstů upínadla" [▶ 187]. Pomocí momentového šroubováku utáhněte šrouby na 3 Nm, jak je popsáno v části "Základní nastavení prstů upínadla pro prsty upínadla FES" [▶ 188].

8.2.5.1 Řešení problémů s robotickým upínacím ramenem s dlouhou osou Z (RGA-Z)

Tab. 33: Tabulka řešení problémů s robotickým upínacím ramenem s dlouhou osou Z (RGA-Z)

Problém/chyba	Možná příčina	Nápravné opatření
Mikrodeska nebyla vyzvednuta	Žádné mikrodesky na nosiči. Prsty upínadla nemohou mikrodesku zvednout.	Položte mikrodesku na nosič. Nastavte polohu upínadla. Vyčistěte prsty upínadla RGA.
	Prsty upínadla jsou kluzké.	Vyčistěte prsty upínadla RGA.
Neobvyklý hluk při pohybu ramena	Díly jsou poškozené nebo opotřebované.	Viz část "Zákaznická podpora" [▶ 199].

8.2.6 Řešení problémů mycího systému

Tab. 34: Tabulka řešení problémů mycího systému

Problém/chyba	Možná příčina	Nápravné opatření
Chybné přetečení a prázdné chyby během běhu procesu	Snímač přetečení není čistý. Snímače nejsou připojeny nebo jsou vadné.	Zkontrolujte, zda je snímač připojen. Viz příručku výrobce snímače.
Pravidelné přetečení mycí stanice	Odpadní čerpadlo je vadné. Trubky odpadního čerpadla jsou zalomené, ucpané nebo poškozené.	Zkontrolujte odpadní potrubí. V případě potřeby vyměňte odpadní trubky. Viz část "Zákaznická podpora" [▶ 199].



Problém/chyba	Možná příčina	Nápravné opatření
Mycí stanicí neprotéká žádná mycí kapalina nebo je mycí systém prázdný	Trubky odpadního čerpadla jsou zalomené, ucpané nebo poškozené.	Zkontrolujte odpadní potrubí. V případě potřeby vyměňte odpadní čerpadlo. Viz část "Zákaznická podpora" [▶ 199].
	Mycí stanice není připojena. Nádoba (nádoby) na mytí je (jsou) prázdná (prázdné) nebo chybí.	Doplňte nebo vyměňte mycí nádobu (nádoby). Správně připojte mycí stanici (MCA). Viz část "Připojení mycí stanice (MCA 384)" [▶ 166].
	Odpadní čerpadlo je vadné.	Zkontrolujte odpadní čerpadlo. V případě potřeby vyměňte odpadní čerpadlo. Viz část "Zákaznická podpora" [▶ 199].
Přetečení mycí stanice	Odpadní trubka je pod hladinou kapaliny v nádobě na odpad.	Použijte mycí nádobu s pevným přívodem mycí trubky.
	Jednorázové hroty nebo řasy blokují mycí stanici.	Vyčistěte mycí stanici. Viz část "Čištění nádoby systémové kapaliny a nádoby na odpad" [▶ 142].
	Odpadní potrubí je zalomené.	Zkontrolujte, zda nejsou trubky zalomené. Viz část "Kontrola potrubí na nádobě systémové kapaliny a nádobě na odpad" [▶ 91].



8.2.7 Řešení problémů s Fluent ID

Tab. 35: Tabulka řešení problémů s Fluent ID

Problém/chyba	Možná příčina	Nápravné opatření
Čárový kód nebyl přečten	Štítek s čárovým kódem nesměřuje ke skeneru.	Vyložte podložku zkumavek, otočte zkumavky tak, aby štítky s čárovým kódem směřovaly doleva. Vložte znovu podložku zkumavek na Fluent.
	Podložka se vkládá příliš rychle.	Vyložte podložku zkumavek a znovu ji pomalu vložte.
	Špatná kvalita štítku.	Zadejte čárový kód ručně nebo nahlaste problém hlavnímu operátorovi.
	Okno skeneru je špinavé.	Vyčistěte okno skeneru. Viz část "Týdenní péče o systém" [▶ 119].
	Reflektor je špinavý.	Vyčistěte reflektor. Viz část "Týdenní péče o systém" [▶ 119].
	Typ čárového kódu nebo délka čárového kódu nejsou pro metodu předem definovány.	Nahlaste problém hlavnímu operátorovi.
Přítomnost zkumavky nebyla zjištěna	Umístění štítku s čárovým kódem příliš nízko na zkumavce.	Nahlaste problém hlavnímu operátorovi.

8.2.8 Řešení problémů se softwarem

Tab. 36: Tabulka řešení problémů se softwarem

Problém/chyba	Možná příčina	Nápravné opatření
Přihlašovací obrazovka uživatele se nezobrazuje, když se to očekává.	Správa uživatelů nebyla ve FluentControl aktivována.	Pro aktivaci správy uživatelů se obraťte na hlavního operátora.
Uživatel se nemůže přihlásit.	Heslo je nesprávné nebo je účet uzamčen.	Pro obnovení hesla nebo účtu se obraťte na hlavního operátora.
Ne všechny servisní úkony byly dokončeny. Upozornění se zobrazí při každém spuštění FluentControl.	Ne všechny předpokládané servisní úkony jsou v konfiguraci přístroje označeny jako dokončené.	Viz část "Zákaznická podpora" [▶ 199].



Problém/chyba	Možná příčina	Nápravné opatření
Dotyková obrazovka nereaguje na dotyk.	Softwarový ovladač není nainstalován.	Pro instalaci ovladačů na instalačním CD a konfiguraci dotykové obrazovky se obraťte na správce počítače.
	Nesprávně nakonfigurované rozhraní dotykové obrazovky.	Otevřete nastavení ovladače dotykové obrazovky a zkontrolujte, zda je dotyková obrazovka správně namapována.
Dotykové rozhraní se nezobrazuje na dotykové obrazovce.	Dotyková obrazovka nebyla při spuštění softwaru zapnutá.	Zapněte přístroj a restartujte software nebo zkontrolujte nastavení dotykového nástroje v konfiguračním systému FluentControl.
Chyba při spuštění FluentControl.	Program FluentControl (SystemSW.exe) je již spuštěn na pozadí (Správce úloh).	Otevřete Správce úloh, zpracujte SystemSW.exe a restartujte FluentControl. Nebo restartujte počítač.
FluentControl nekomunikuje s připojenými hardwarovými zařízeními.	FluentControl není správně nakonfigurován pro komunikaci s hardwarovými zařízeními.	Obraťte se na osobu odpovědnou za konfiguraci systému, aby aktivovala stav I/ O hardwarových zařízení.



8.3 Činnosti při řešení problémů

8.3.1 Připojení mycí stanice (MCA 384)

Chcete-li připojit mycí stanici ramena MCA 384, postupujte následovně:

- 1. Vypněte přístroj.
- 2. Připojte mycí trubku (A) a odpadní trubku (B) k mycí stanici.



3. Připojte mycí trubku (A) a odpadní trubku (B).







4. Připojte všechny trubky (C) k řídicí jednotce mycí stanice ramena MCA 384.

5. Obraťte se na hlavního operátora a zkontrolujte připravenost k provozu.

8.3.2 Pozice laboratorního vybavení

Chcete-li zajistit správné umístění laboratorního vybavení v soupravě pro přesný přístup k ramenu, postupujte následovně:

1. Umístěte laboratorní vybavení (A) na soupravu (B).







2. Jemně přitlačte laboratorní vybavení ke statickému polohovacímu zařízení (C).

- 3. Posuňte diagonální posuvné polohovadlo (E) směrem k laboratornímu vybavení nebo od něj, aby se přesně přizpůsobil laboratornímu vybavení.
- 4. Posuňte vertikální a horizontální posuvné polohovadlo (D, F) směrem k laboratornímu v vybavení nebo od něj, abyste laboratorní vybavení upevnili.
- 5. Zvedněte laboratorní vybavení z hnízda.

Dbejte na to, aby při nasazování nebo snímání desky nedocházelo ke tření.

8.3.3 Kontrola řadového filtru (FCA se vzduchem)

Na každém kanálu je nainstalován řídicí systém, který chrání pipetovací kanály vzduchového ramena FCA před nadměrnou aspirací kapaliny.

- Rutinní údržba FCA se vzduchem zahrnuje kontrolu řadových filtrů, která odhalí mokré, poškozené nebo špatně umístěné filtry a chybějící filtry.
- Spusťte metodu Rutinní údržba FCA se vzduchem za účelem kontroly řadového filtru uvnitř kuželu DiTi pipetovacího kanálu vzduchového ramena FCA.

V případě chyby je potřeba vyměnit řadový filtr. Viz část "Výměna řadového filtru (FCA se vzduchem)" [▶ 168].

8.3.4 Výměna řadového filtru (FCA se vzduchem)

Chcete-li vyměnit řadový filtr, postupujte následovně:

- ✓ Dekontaminovaný kužel s jednorázovým hrotem.
- ✓ Kužel DiTi byl odebrán z kanálu. Viz část "Demontáž kuželu DiTi (FCA se vzduchem)" [▶ 170].





- 1. Propíchněte řadový filtr (B) do strany nástrojem pro demontáž filtru (C).
- Vypačte řadový filtr pomocí nástroje na demontáž filtru. Povšimněte si, že filtr může být znečištěn procesními kapalinami.
- 3. Zlikvidujte řadový filtr.



- Vyčistěte kužel DiTi (A) alkoholem. Kužel DiTi musí být před opětovnou montáží suchý.
- 5. Umístěte nový řadový filtr na čistý a rovný povrch.
- 6. Vtlačte řadový filtr do kuželu DiTi.

Řadový filtr nesmí vyčnívat z kuželu DiTi.

7. Zkontrolujte řadový filtr podle metody definované vaším hlavním operátorem.



8.3.5 Demontáž kuželu DiTi (FCA se vzduchem)

Chcete-li vyjmout kužel DiTi (FCA se vzduchem), postupujte takto:

- ✓ K dispozici je klíč na kužel DiTi.
- 1. Vypněte přístroj.
- 2. Otevřete přední bezpečnostní panel.
- 3. Ručně zvedněte všechny tyče Z do nejvyšší polohy.
- 4. Přesuňte všechny tyče Z směrem k přední části přístroje.
- 5. Roztáhněte tyče Z co nejvíce do šířky.
- 6. Přidržte adaptér hrotu (D) a trubku pro vyhazování hrotu (C).
- 7. Odšroubujte kužel DiTi (A) pomocí klíče na kužel DiTi (B).



8. Opatrně stáhněte kužel DiTi.



V některých případech může být ke kuželu DiTi (A) stále připojena vyhazovací trubice hrotu (C) nebo válec adaptéru (B). Viz část "Montáž vyhazovací trubice DiTi (FCA se vzduchem)" [▶ 171].



8.3.6 Montáž vyhazovací trubice DiTi (FCA se vzduchem)

Chcete-li namontovat vyhazovací hadici DiTi (Air FCA), postupujte takto:

- ✓ Vyhazovací hadice DiTi byla vyjmuta podle pokynů.
- ✓ K dispozici je klíč na kužel DiTi.
- 1. Vložte těsnicí pouzdro (B) do válce adaptéru (A).









2. Ke smontovanému válci přišroubujte kužel DiTi (C). Zkontrolujte, že černý Okroužek není vidět, jak je znázorněno na obrázku dole.

3. Zasuňte vyhazovací hadici špiček (D) stranou bez zářezu do sestaveného válce, jak je znázorněno níže.



8.3.7 Montáž kuželu DiTi (FCA se vzduchem)

Při montáži kužele DiTi vzduchového ramena FCA postupujte následovně:

- ✓ Kužel DiTi je kompletně sestaven: Viz část "Montáž vyhazovací trubice DiTi (FCA se vzduchem)" [▶ 171].
- ✓ K dispozici je klíč na kužel DiTi.
- 1. Zasuňte válec adaptéru do trubky vyhazovače hrotů (C).
- 2. Přidržte adaptér hrotu (D) a trubku pro vyhazování hrotu (C).





3. Našroubujte kužel DiTi (A) pomocí klíče na kužel DiTi (B).

4. Spusťte metodu Rutinní údržba FCA se vzduchem .



8.3.8 Demontáž doplňku DiTi (FCA)

Chcete-li odstranit doplněk DiTi, postupujte takto:

- ✓ Klíč na kužel DiTi
- 1. Vypněte přístroj.
- 2. Otevřete přední bezpečnostní panel.
- 3. Ručně zvedněte všechny tyče Z do nejvyšší polohy.
- 4. Přesuňte všechny tyče Z směrem k přední části přístroje.
- 5. Roztáhněte tyče Z co nejvíce do šířky.
- 6. Přidržte adaptér hrotu (D) a trubku pro vyhazování hrotu (C).
- 7. Odšroubujte kužel DiTi (A) pomocí klíče na kužel DiTi (B).



8. Opatrně stáhněte kužel DiTi.

8.3.9 Montáž doplňku DiTi (FCA)

Chcete-li namontovat doplněk DiTi, postupujte takto:

- 1. Našroubujte válec adaptéru (B) do trubky vyhazovače hrotů (C).
- 2. K adaptéru válci přišroubujte kužel DiTi (A). Použijte klíč na kužel DiTi.







3. Zatlačte trubku na plastovou jehlu, dokud se trubka pevně nepřipojí k doplňku DiTi.

- 4. Přidržte adaptér hrotu (D) a trubku pro vyhazování hrotu (C).
- 5. Našroubujte kužel DiTi (A) pomocí klíče na kužel DiTi (B).



8.3.10 Kontrola fixních hrotů

Při kontrole fixních hrotů postupujte následovně:

OZNÁMENÍ

Nepřesnost pipetování a chyby při detekci kapalin!

Ohnutý nebo poškozený povlak hrotů způsobují nepřesnost pipetování a chyby při detekci kapalin.

• Nikdy nepracujte s poškozenými nebo ohnutými hroty.



- 1. Vypněte přístroj.
- 2. Otevřete přední bezpečnostní panel.
- 3. Zkontrolujte fixní hroty.
- 4. Zkontrolujte povlak fixního hrotu zrcátkem.

Ujistěte se, že fixní hroty nejsou ohnuté. Pokud je povlak fixního hrotu poškozen nebo je fixní hrot ohnutý, je potřeba jej vyměnit. Viz část "Demontáž fixních hrotů" [> 176].

8.3.11 Demontáž fixních hrotů

Při demontáži fixních hrotů postupujte následovně:

- ✓ Fixní hroty byly vyčištěny. Viz část "Tabulky péče o systém" [▶ 114].
- \checkmark Fixní hroty byly zkontrolovány. Viz část "Kontrola fixních hrotů" [> 175].



- С Pojistná matice
- D
- Е Pipetovací potrubí
- 1. Vypněte přístroj.
- 2. Otevřete přední bezpečnostní panel.
- 3. Ručně zvedněte všechny tyče Z (A) do nejvyšší polohy.
- 4. Roztáhněte tyče Z co nejvíce do šířky.
- Pokud je nainstalován nastavitelný fixní hrot, povolte čtyři šrouby pro nastavení 5. hrotu.
- 6. Odšroubujte pojistnou matici (C) a druhou rukou přidržujte fixní hrot těsně pod pojistnou maticí.
- 7. Odstraňte pojistnou matici (C) a posuňte ji podél osy hrotu. Zabraňte kontaktu pojistné matice s povlakem hrotu.
- 8. Pokud je hrot (D) nastavitelný, otočte pojistnou matici (C) vzhůru nohama na čistý povrch a vyjměte O-kroužek a podložku.



- Pokud je kanál vybaven doplňkem s malým objemem, odšroubujte přírubu na horní straně elektromagnetického ventilu, abyste uvolnili pipetovací trubku (E) procházející tyčí Z (A).
- 10. Vytáhněte pipetovací trubku (E) o určitou vzdálenost (a) z adaptéru hrotu (B) tahem za hrot (D).

Použijte suché smirkové plátno pro lepší uchopení pipetovací trubky – nikoli hrotu.

8.3.12 Montáž fixních hrotů



Při montáži fixních hrotů postupujte následovně:

1. Opatrně vytáhněte pipetovací trubku o délku přibližně 25 mm (1 palec) (a) z adaptéru hrotu.

Použijte malý kousek smirkového plátna, abyste trubku na konci lépe uchopili. Pokud byl hrot nainstalován již dříve, odřízněte z pipetovací trubky přibližně 5 mm (0,2 palce) (b) pomocí ostrého nože, abyste dosáhli rovného řezu.





- Nastavitelná pojistná matice K Nastavovací šroub hrotu
- 2. V případě hrotů Te-PS nebo hrotů s malým objemem:

Pomocí rozšiřovače trubek Te-PS (G) rozšiřte konec trubky tak, že rozšiřovač trubek Te-PS zatlačíte až po rukojeť do trubky a zároveň otáčíte nástrojem. Zatímco je trubka stále široká, zatlačte hrot Te-PS do konce trubky přibližně o 4 mm (0,16 palce).

3. Nasaďte pojistnou matici na hrot.

J

Pokud je hrot nastavitelný (např. Te-PS), nasaďte pojistnou matici na podložku (H) a O-kroužek (I).

OZNÁMENÍ Vyvarujte se kontaktu s jemným koncem hrotu a jeho povlakem.

- 4. Vložte hrot a pipetovací trubku do adaptéru hrotu.
- 5. Našroubujte pojistnou matici na adaptér hrotu a utáhněte ji.

Pokud je hrot nastavitelný (např. Te-PS), utáhněte pojistnou matici tak, aby čtyři šrouby pro nastavení hrotu (K) svíraly úhel 45° se souřadnicovým systémem plošiny X/Y.

- 6. Vyčistěte fixní hroty. Viz část "Konec dne" [▶ 116].
- 7. Proveďte test přesnosti pipetování podle definice hlavního operátora.





8.3.13 Demontáž propichovacích hrotů

Při demontáži propichovacího hrotu postupujte následovně:

- ✓ Přístroj je vypnutý.
- 1. Otevřete přední bezpečnostní panel.
- 2. Ručně zvedněte všechny tyče Z do nejvyšší polohy.
- 3. Přesuňte všechny tyče Z směrem k přední části přístroje.
- 4. Roztáhněte tyče Z co nejvíce do šířky.





5. Zakryjte propichovací hroty ochrannými pomůckami. Začněte nejzadnějším propichovacím hrotem.

6. Odšroubujte pojistnou matici, druhou rukou přidržujte propichovací hrot těsně pod pojistnou maticí.




7. Tažením za hrot vytáhněte pipetovací trubku asi 25 mm z adaptéru hrotu. Při vytahování držte propichovací hrot na jeho horním konci.

- 8. Vytáhněte propichovací hrot z trubky a druhou rukou trubku přidržujte.
- 9. Neodstraňujte ochranu propichovacího hrotu. Vyhoďte ji i s propichovacím hrotem do kontejneru na biologický odpad.







8.3.14 Montáž propichovacích hrotů

- Přístroj je vypnutý.
- ✓ Hlavní operátor je dostupný.
- 1. Otevřete přední bezpečnostní panel.
- 2. Ručně zvedněte všechny tyče Z do nejvyšší polohy.
- 3. Přesuňte všechny tyče Z směrem k přední části přístroje.
- 4. Roztáhněte tyče Z co nejvíce do šířky.



5. Otevřete obal propichovacího hrotu. Ochranu hrotu (F) neodstraňujte.



Pořadí instalace propichovacích hrotů: zezadu dopředu

- 6. Opatrně vytáhněte pipetovací trubku o délku přibližně 25 mm z adaptéru hrotu.
- 7. Zasuňte prázdný kuželový konec propichovacího hrotu do konce trubky.



8. Vložte propichovací hrot a pipetovací trubku do adaptéru hrotu.





9. Našroubujte pojistnou matici na adaptér hrotu a utáhněte ji rukou.

- 10. Mírně otevřete pojistnou matici. Posuňte ochranu hrotu mírně dolů, abyste získali přístup k dříku propichovacího hrotu. Ochranu hrotu ještě zcela neodstraňujte.
- 11. Otočte propichovací hrot tak, aby otvor hrotu směřoval k přední straně přístroje. Jednou rukou udržujte propichovací hrot v této orientaci a druhou rukou dotáhněte pojistnou matici.
- 12. Zkontrolujte, zda všechny otvory hrotu směřují k přední straně přístroje.







13. Po instalaci všech propichovacích hrotů odstraňte všechny ochrany hrotů. Začněte nejzadnějším propichovacím hrotem.

- 14. Pro vynulování čítače v FluentControl se obraťte na hlavního operátora.
- 15. Kontaktujte hlavního operátora pro provedení testu soupravy QC. Viz Referenční dokumenty.
- 16. Spusťte metodu Únik propichování FCA .
- 17. Proveďte test přesnosti pipetování (doporučení: použijte sadu QC) podle definice hlavního operátora.



8.3.15 Vytahování zaseknutých propichovacích hrotů



V případě, že se propichovací hrot zasekne tak, že jej nelze vytáhnout pomocí softwarových příkazů, je nutné jej vyjmout ručně.

Při vytahování propichovacích hrotů postupujte následovně:

- ✓ Přístroj je vypnutý.
- 1. Otevřete přední bezpečnostní panel.
- 2. Ručně zvedněte všechny zasunuté tyče Z do nejvyšší polohy.
- Zakryjte všechny zasunuté propichovací hroty ochrannými kryty propichovacích hrotů (A). Začněte nejzadnějším propichovacím hrotem.





4. Umístěte nástroj pro odstranění hrotu vedle zaseknutého hrotu na pevný a stabilní povrch a nasaďte jej pod pojistnou matici.



- 5. Otáčejte knoflíkem nástroje pro odstranění propichovacího hrotu, dokud není hrot zcela vytažen.
- 6. Otočte knoflíkem v opačném směru a snižte retraktor asi o 1 cm.
- 7. Vyjměte nástroj pro odstranění propichovacího hrotu.
- Zaseknutý hrot je nyní vytažen.8. Vyčistěte nástroj pro odstranění propichovacího hrotu alkoholem.
- 9. Zkontrolujte, zda není propichovací hrot poškozen (např. ohnutý propichovací hrot, poškozený hrot).
- 10. Pokud je propichovací hrot poškozený, vyměňte jej. Viz část "Demontáž propichovacích hrotů" [▶ 179] a část "Montáž propichovacích hrotů" [▶ 182].
- 11. Jednou rukou držte pojistnou matici a druhou rukou sejměte všechny ochrany propichovacího hrotu. Začněte nejzadnějším propichovacím hrotem.
- 12. Vyčistěte nástroj pro odstranění propichovacího hrotu alkoholem.

Also see about this

- "Demontáž propichovacích hrotů" > 179
- "Montáž propichovacích hrotů" > 182

8.3.16 Kontrola vyrovnání prstů upínadla

Po srážce nebo při nasazení náhradních prstů upínadla může být nutné provést seřízení prstů upínadla. To platí pro všechna ramena používající upínadla s prsty upínadel.



Chybné vyrovnání po srážce:

- Analyzujte situaci.
- Vyhodnoťte možné příčiny srážky nebo příčinu špatného nastavení prstu, například špatně nastavenou zásuvku čtečky, podložku, nesprávně naučený/ umístěný hotel nebo jiný segment.
- Vyberte níže uvedený postup na základě požadavků na přesnost.



- Pokud prsty upínadla nepotřebují splňovat nadprůměrnou přesnost, proveďte základní vyrovnání. Viz část "Základní nastavení prstů upínadla pro prsty upínadla FES" [▶ 188], nebo část "Základní nastavení prstů upínadla pro fixní prsty upínadla" [▶ 189].
- Pokud musí prsty upínadla splňovat pokročilé požadavky (odchylka Z < ±0,2 mm), proveďte postup pokročilého vyrovnání prstů upínadla. Viz část "Pokročilé nastavení prstů upínadla pro prsty upínadla FES" [▶ 190], nebo "Pokročilé nastavení prstů upínadla pro fixní prsty upínadla" [▶ 190].
- 3. Prsty upínadla lze namontovat pomocí dvou různých typů šroubů: a) šroub M4 × 12 s šesticípou hvězdicí, utažený momentem 3 Nm.



b) inbusový šroub M4 × 12 v kombinaci s napínací podložkou (dodržujte polohu podle obrázku níže), utažený momentem 3,5 Nm.





Pokud nemáte k dispozici momentový šroubovák, utahujte šroub tak dlouho, dokud nebude podložka stlačena naplocho a odpor se nezvýší. Poté dotáhněte o ¹/₁₂ otáčky. To odpovídá přibližně 3,5 Nm.

8.3.17 Základní nastavení prstů upínadla pro prsty upínadla FES

To platí pro všechna ramena používající upínadla s prsty upínadel.

Při základním vyrovnání postupujte následovně:

- ✓ Chybné vyrovnání je jasně viditelné.
- Není vyžadována nadprůměrná přesnost.
- ✓ K dispozici je momentový šroubovák (s možností nastavení 3 nebo 3,5 Nm). Pokud není k dispozici momentový šroubovák:
 Šroub s šesticípou hvězdicí: pevně utáhněte šrouby, ale nepoužívejte nadměrnou sílu.
 Šroub s vnitřním šestihranem: viz část "Kontrola vyrovnání prstů upínadla"
 [▶ 187].



- 1. Odstraňte prst upínadla z hlavy upínadla.
- 2. Povolte šroub mezi prstem upínadla a adaptérem prstu FES.
- Přitlačte prst upínadla k hornímu a zadnímu dorazu adaptéru, jak je znázorněno na obrázku níže, a utáhněte šroub momentovým šroubovákem (3 nebo 3,5 Nm).



8.3.18 Základní nastavení prstů upínadla pro fixní prsty upínadla

Při základním vyrovnání postupujte následovně:

- ✓ Chybné vyrovnání je jasně viditelné.
- ✓ Není vyžadována nadprůměrná přesnost.
- ✓ K dispozici je momentový šroubovák (s možností nastavení 3 nebo 3,5 Nm).
- 1. Povolte upevňovací šroub (A).





2. Přitlačte prst upínadla k hornímu a zadnímu dorazu montážního prvku hlavy upínadla a utáhněte šroub momentovým šroubovákem (3 nebo 3,5 Nm).

8.3.19 Pokročilé nastavení prstů upínadla pro prsty upínadla FES

To platí pro všechna ramena používající upínadla s prsty upínadel.

- ✓ K dispozici je momentový šroubovák (s možností nastavení 3 nebo 3,5 Nm). Pokud není k dispozici momentový šroubovák:
 Šroub s šesticípou hvězdicí: pevně utáhněte šrouby, ale nepoužívejte nadměrnou sílu.
 Šroub s vnitřním šestihranem: viz část "Kontrola vyrovnání prstů upínadla"
 [▶ 187].
- 1. Pomocí nástroje přemístění přesuňte výšku Z do výšky přibližně 3 mm nad pracovní stůl.
- 2. Povolte šroub mezi prstem upínadla a adaptérem prstu FES.



- 3. Ujistěte se, že je adaptér prstů FES pevně připojen k hlavě upínadla. *Adaptéry prstů jsou v jednom směru přidržovány magnetem.*
- Pomocí nástroje pro přesouvání přesuňte výšku Z do výšky 0 mm nad pracovní stůl.
 Poslední desetinu milimetru přesouvejte pomalu.
 Poznámka: Pokud nevíte, jak přistupovat k nástroji přemístění nebo jak jej ovládat, obraťte se na hlavního operátora.
- 5. Přitiskněte prst upínadla k hlavě upínadla a k referenčnímu povrchu, jak je znázorněno na obrázku, a utáhněte šroub utahovacím momentem 3 nebo 3,5 Nm.
- Nastavení zkontrolujte ručním otočením hlavy o 90°, 180°, 270°. Chybné vyrovnání v různých polohách znamená nesprávné nastavení hlavy nebo ramena. V tomto případě je třeba, aby vyrovnání zkontroloval technik FSE.

8.3.20 Pokročilé nastavení prstů upínadla pro fixní prsty upínadla

✓ K dispozici je momentový šroubovák (s možností nastavení 3 nebo 3,5 Nm).



- 1. Pomocí nástroje přemístění přesuňte výšku Z do výšky přibližně 3 mm.
- 2. Povolte šroub mezi prstem upínadla a hlavou upínadla.
- Pomocí nástroje přemístění přesuňte výšku Z do výšky 0 mm. Poslední desetinu milimetru se pohybujte pomalu.
 Poznámka: Pokud nevíte, jak přistupovat k nástroji přemístění nebo jak jej ovládat, obraťte se na hlavního operátora.
- 4. Přitiskněte prst upínadla k hlavě upínadla a k referenčnímu povrchu a utáhněte šroub utahovacím momentem 3 nebo 3,5 Nm.
- 5. Nastavení zkontrolujte ručním otočením hlavy o 90°, 180°, 270°. Chybné vyrovnání v různých polohách znamená nesprávné nastavení hlavy nebo ramena. V tomto případě je třeba, aby vyrovnání zkontroloval technik FSE.



8.3.21 Kontrola segmentu

Zkontrolujte, zda je segment uzavřen.



Obr. 69: Segment je uzavřený



Obr. 70: Segment je otevřený



8.3.22 Demontáž polohovacích kolíků

Při demontáži polohovacích kolíků postupujte následovně:

OZNÁMENÍ

Havárie nebo chyba procesu!

Havárie a další chyby procesu mohou být důsledkem nepřesného umístění prvků na segmentu plošiny v důsledku uvolněných polohovacích kolíků.

- Produkt Fluent nepoužívejte, pokud chybí polohovací kolíky.
- ✓ Polohovací kolíky jsou zlomené.
- 1. Nasaďte odstraňovač kolíků na polohovací kolík.







2. Zvedněte rukojeť odstraňovače kolíků a vytáhněte kolík ze segmentu plošiny.

8.3.23 Výměna blokovacích kolíků a polohovacích kolíků

OZNÁMENÍ

Havárie nebo chyba procesu!

Havárie a další chyby procesu mohou být důsledkem nepřesného umístění prvků na segmentu plošiny v důsledku uvolněných polohovacích kolíků.

- Produkt Fluent nepoužívejte, pokud chybí polohovací kolíky.
- ✓ K dispozici jsou polohovací kolíky uvedené v tabulce péče o systém.
- Odstraňte segment z plošiny: Viz část "Odebírání segmentů" [▶ 83].
- 2. Zatlačte nový blokovací kolík (B) do otvoru (A).



- 3. Zatlačte nový polohovací kolík (C) do otvoru (A).

Umístěte segment na plošinu:
 Viz část "Umístění segmentů" [▶ 82].



9 Balení, vybalování, přeprava, skladování a likvidace

Tato kapitola uvádí regulační informace o recyklaci a etiketách na obalech, které je potřeba dodržovat.

OZNÁMENÍ

Zabraňte poškození nekvalifikovaným a neoprávněným personálem!

Balení, vybalování, přepravu a skladování smí provádět pouze pracovníci společnosti Tecan nebo pracovníci pověření společností Tecan!
Viz "Zákaznická podpora" [> 199].

Informace o přemístění přístroje naleznete v části "Přemístění přístroje na skříni v rámci laboratoře" [▶ 124].

9.1 Etikety na obalu

Správné a úplné označení obalů pomáhá předcházet nesprávné manipulaci, nehodám, nesprávnému dodání, ztrátě hmotnosti a poškození během skladování.

Symbol	Význam	Popis
	Recyklujte	Obalový materiál lze recyklovat. Nelikvidujte jako domácí odpad. Informace o materiálu použitém pro tento obal jsou uvedeny pod symbolem.
<u>11</u>	Touto stranou nahoru	Dbejte na to, aby byl obal přepravován a skladován horní stranou, označenou šipkami, nahoře. Nepřevracejte.
	Udržujte v suchu	Zajistěte, aby se obal během přepravy a skladování nenamočil.
	Křehké	S balením zacházejte opatrně. Uvnitř je křehké zboží.
*	Chraňte před slunečním zářením	Zajistěte, aby obal nebyl během přepravy a skladování vystaven teplu. Chraňte před silným slunečním zářením.
	Neukládejte na sebe	Balení neukládejte na sebe. Obal není navržen tak, aby unesl dodatečnou hmotnost.

Tab. 37: Symboly na obalu



9.2 Likvidace

Tato část uvádí informace o předpisech týkajících se recyklace, které je třeba dodržovat.

OZNÁMENÍ

Recyklujte v souladu s platnými právními předpisy!

Dodržujte zákony platné ve vaší zemi pro recyklaci.

9.2.1 Místní požadavky Evropské unie

Evropská komise vydala směrnici o odpadních elektrických a elektronických zařízeních (OEEZ; 2012/19/EU).

Od srpna 2005 jsou výrobci zodpovědní za zpětný odběr a recyklaci elektrických a elektronických zařízení.

Označení	Vysvětlení
	Negativní dopady na životní prostředí spojené se zpracováním odpadu.
	 S elektrickými a elektronickými zařízeními nezacházejte jako s netříděným komunálním odpadem.
	 Odpadní elektrická a elektronická zařízení sbírejte odděleně.

9.2.2 Místní požadavky Čínské lidové republiky

Označení pro omezení používání nebezpečných látek v elektronických a elektrických výrobcích

Norma pro elektronický průmysl Čínské lidové republiky SJ/T11364-2014 Označení pro omezení používání nebezpečných látek v elektronických a elektrických výrobcích vyžaduje označení pro omezení používání nebezpečných látek v elektronických a elektrických výrobcích.

V souladu s požadavky uvedenými v SJ/T11364-2014 jsou všechny elektronické a elektrické výrobky Tecan prodávané v Čínské lidové republice označeny značkou pro omezení používání nebezpečných látek.

Označení	Vysvětlení
25	Toto označení znamená, že tento elektronický výrobek obsahuje určité nebezpečné látky a může být bezpečně používán po dobu používání šetrného k životnímu prostředí, ale po uplynutí doby používání šetrného k životnímu prostředí musí být zařazen do recyklačního systému.



9.2.3 Ostatní požadavky

Označení	Vysvětlení
Hg	 Tato lampa obsahuje rtuť Recyklujte nebo likvidujte podle platných místních zákonů.



10 Zákaznická podpora

Tato kapitola vysvětluje, které soubory a informace společnost Tecan vyžaduje k prvnímu posouzení problému.

10.1 Kontakty

Obraťte se na místního distributora či dovozce nebo na jednu z níže uvedených adres.

Podívejte se také na naši domovskou stránku na adrese: www.tecan.com

Tab. 38: Kontakty na zákaznickou podporu

Země/region	Adresa	Telefon/fax/e-	mail
Asie	Tecan Asia Pte Ltd. 18 Boon Lay Way, #10-106 TradeHub 21 Singapore 609966 Singapore	Telefon Fax E-mail	+65 6444 1886 +65 6444 1836 tecan@tecan.com.sg
Austrálie Nový Zéland Tichomořské ostrovy	Tecan Australia Pty Ltd Unit 2, 475 Blackburn Road Mount Waverly VIC 3149 Austrálie	Telefon Telefon Fax E-mail	Bezplatná telefonní linka: 1300 808 403 +61 3 9647 4100 +61 3 9647 4199 helpdesk-aus@tecan.com
Rakousko	Tecan Sales Austria GmbH Untersbergstrasse 1a 5082 Grödig Rakousko	Telefon Fax E-mail	+43 6246 8933 256 +43 6246 72770 helpdesk-at@tecan.com
Belgie	Tecan Benelux B.V.B.A. Mechelen Campus Schaliënhoevedreef 20A 2800 Mechelen Belgie	Telefon Fax E-mail	+32 15 42 13 19 +32 15 42 16 12 tecan-be@tecan.com
Čína	Tecan (Shanghai) Laboratory Equipment Co., Ltd. 1F, T 15-4, #999, Ningqiao Road, Pilot Free Trade Zone, Shanghai, PRC, 201206	Telefon E-mail	+86 40 0821 38 88 helpdesk-cn@tecan.com



Země/region	Adresa	Telefon/fax/e-	mail
Dánsko	Tecan Denmark, Filial af Tecan Nordic AB, Sverige Lejrvej 29 3500 Værløse Dánsko	Telefon E-mail	+46 8 7503940 info-dk@tecan.com
Francie	Tecan France S.A.S.U Tour Swiss Life 1 bd Marius Vivier Merle F- 69 003 Lyon Francie	Telefon Fax E-mail	+33 4 72 76 04 80 +33 4 72 76 04 99 helpdesk-fr@tecan.com
Německo	Tecan Deutschland GmbH Werner-von-Siemens-Straße 23 74564 Crailsheim Německo	Telefon Fax E-mail	+49 1805 8322 633 nebo +49 1805 TECAN DE +49 7951 9417 92 helpdesk-de@tecan.com
Itálie	Tecan Italia, S.r.I. Via Brescia, 39 20063 Cernusco Sul Naviglio (MI) Itálie	Telefon Fax E-mail	+39 800 11 22 91 +39 (02) 92 72 90 47 helpdesk-it@tecan.com
Japonsko	Tecan Japan Co., Ltd. Kawasaki Tech Center 580-16, Horikawa-cho, Saiwai-ku Kawasaki, Kanagawa 212-0013 Japonsko	Telefon Fax Telefon E-mail	+81 44 556 7311 (Kawasaki) +81 44 556 7312 (Kawasaki) +81(0) 6305 8511 (Osaka) helpdesk-jp@tecan.com
Nizozemsko	Tecan Benelux B.V.B.A. Industrieweg 30 NL-4283 GZ Giessen Nizozemsko	Telefon Fax E-mail	+31 20 708 4773 +31 183 44 80 67 helpdesk.benelux @tecan.com
Skandinávie	Tecan Nordic AB Sveavägen 159, 1tr SE-113 46 Stockholm Švédsko	Telefon Fax E-mail	+46 8 750 39 40 +46 8 750 39 56 info@tecan.se
Jižní Korea	Tecan Korea Ltd. 149 Gasan digital 1-ro Geumcheon-gu Seoul Jižní Korea	Telefon E-mail	+82-2-818-3301 helpdesk-kr@tecan.com



Země/region	Adresa	Telefon/fax/e-	mail
Španělsko Portugalsko	Tecan Ibérica Instrumentación S.L. C/ Lepanto 151 Bajos E-08013 Barcelona Španělsko	Telefon E-mail	+34 93 595 25 31 helpdesk-sp@tecan.com
Švýcarsko	Tecan Schweiz AG Seestrasse 103 8708 Männedorf Švýcarsko	Telefon Fax E-mail	+41 44 922 82 82 +41 44 922 89 23 helpdesk-ch@tecan.com
Spojené království	Tecan UK Ltd. Theale Court 11-13 High Street Theale, Reading, RG7 5AH Spojené království	Telefon Fax E-mail	+44 118 930 0300 +44 118 930 5671 helpdesk-uk@tecan.com
USA	Tecan US, Inc. 9401 Globe Center Drive, Suite 140, Morrisville, NC 27560 USA	Telefon Fax Telefon E-mail	+1 919 361 5200 +1 919 361 5201 Bezplatná telefonní linka v USA: +1 800 TECAN US nebo +1 800 832 2687 helpdesk-us@tecan.com
USA (Tecan Systems)	Tecan Systems, Inc. 2450 Zanker Road San Jose, CA 95131 USA	Telefon Fax E-mail	+1 408 953 3100 Bezplatná telefonní linka: +1 800 231 0711 +1 408 953 3101 helpdesk-sy@tecan.com



Zkratky

ADT

Air displacement technology – technologie vytlačování vzduchu

ASM

Application Software Manual – Příručka aplikačního softwaru

CE

Conformité Européenne

cLLD

Capacitive Liquid Level Detection – kapacitní detekce hladiny kapaliny

CNS

Common Notification System – Společný oznamovací systém

DiTi

DiTi – jednorázová špička

EMC

Electromagnetic Compatibility – elektromagnetická kompatibilita

EN

Evropská norma

FCA

Flexible Channel Arm – flexibilní rameno kanálu

FCA s kapalinou

Flexibilní rameno kanálu s kapalinovým systémem

FCA se vzduchem

Flexibilní rameno kanálu se vzduchovým systémem

FES

Systém výměny prstů

FSE

Servisní inženýr v terénu

GLP

Správná laboratorní praxe

HEPA

Vysoce účinný filtr pevných částic

IEC

International Electrotechnical Commission – Mezinárodní elektrotechnická komise

IQ

Kvalifikace pro instalaci

ISO

International Organization for Standardization – Mezinárodní organizace pro normalizaci

LED

Light emitting diode – světelná dioda

MCA

Vícekanálové rameno

MCA

Vícekanálové rameno

MET

Registrovaná značka Eurofins EE jako celostátně uznávané zkušební laboratoře

MCH

Multiple channel head – vícekanálová hlava

MIO

Doplněk monitorovaných inkubátorů

•TECAN•

MP

Mikrodeska

NRTL

Celostátně uznávaná zkušební laboratoř

OM

Operating Manual – návod k obsluze

OQ

Provozní kvalifikace

PC

Personal computer - osobní počítač

PP

Polypropylen

rcf

relativní odstředivá síla

RF

Radio Frequency - rádiová frekvence

RGA

Robotic Gripper Arm – robotické upínací rameno

RGA long Z

Robotic Gripper Arm long height – robotické upínací rameno s dlouhou výškou

RGA standard Z

Robotic Gripper Arm standard height – robotické upínací rameno se standardní výškou

RUO

Jen pro výzkumné použití

RWP

RapidWash Pump – čerpadlo RapidWash

SN

Sériové číslo

Te-Shake

Třepačka Tecan

Te-VacS

Vakuový separátor Tecan

USB

Universal Serial Bus – univerzální sériová sběrnice

WEEE

Waste Electrical and Electronic Equipment – odpadní elektrická a elektronická zařízení

WRC

Mycí a doplňovací centrum