

# **Operating Manual**

# Fluent®



Titel:	Bedieningshandleiding Fluent			Onderdeel- nummer:	30199936.03
ID:	399706, nl, V2.7			Vertaald uit het:	n.v.t.
Versie:	Revisie:	Uitgave:	Documentgesch	iedenis:	·
1	0	30-08-2017	Eerste editie		
1	1	18-09-2017	2.3 Toepassings	srisico's	
1	2	07-11-2017	1.3 Beoogd geb	oruik	
1	3	05-03-2018	1.3 Beoogd geb	oruik	
1	4	16-04-2018	3.3 Afmetingen	en gewichten	
1	5	20-07-2018	4.6.2 FCA-grijp	er	
1	6	26-02-2019	2.2 Algemene v	eiligheidsinformat	ie
1	7	22-10-2020	Informatie over Mix & Pierce toegevoegd in ragraaf 1.8.3, 2.5, 2.8.1, 4.6, 4.6.3, 6.6, 7 7.4.2, 7.5.4, 8.2.3, 8.3.13–15. Informatie of buisrotator toegevoegd in paragraaf 4.6.3, 9 7.4.1, 7.4.2, 7.5.5, 7.5.6, 8.2.3, 8.3.14. P graaf 4.6.4 en 7.5.24 toegevoegd. Paragras 2.5, 2.9, 2.10, 4.2.2, 4.6, 6.3.5, 7.1, 7.4, 7.5.16, 8.2.4 en 8.3.13–20 bijgewerkt.		gevoegd in pa- 6.3, 6.6, 7.4.1, nformatie over aaf 4.6.3, 6.5.3, , 8.3.14. Para- d. Paragraaf , 7.1, 7.4, werkt.
1	8	01-12-2020	Paragraaf 2.2 e	n 2.8 bijgewerkt.	
1	9	15-02-2021	1 Paragraaf 2.7.2, 7.1, 7.2, 7.4.2, 7.5.15, 7 en 8.2.3 bijgewerkt, paragraaf 2.7.5, 4.6.4 7.5.25 toegevoegd.		7.5.15, 7.5.24 7.5, 4.6.4 en
1	10	25-11-2021	Paragraaf 3.7, 6.5.1 bijgewerkt.		
2	0	02-03-2022	Paragraaf 3.1 b	ijgewerkt.	
2	1	14-04-2022	Paragraaf 5.2.5 6.5.1 bijgewerkt	, 6.6 toegevoegd	, paragraaf
2	2	16-11-2022	Paragraaf 4.6.5 3.4 bijgewerkt.	toegevoegd, par	agraaf 2.2 en
2	3	15-03-2023	Paragraaf 1.8.2	, 2.4, 3.4 en 7.4.	.2 bijgewerkt.
2	4	30-03-2023	Paragraaf 2.4, 2 4.3.5.3, 7.4.2 e	2.8.1.1, 3.3, 4.3. n 7.4.4 bijgewerk	2, 4.3.3, 4.3.4, kt.

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

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

2	5	16-01-2024	Paragraaf 1.9, 2.2, 2.7.1, 2.8.1, 6.2, 7, 7.2.1, 7.2.2, 7.4.2, 7.5.14, 8.2.2 en 8.3.6 bijgewerkt.
2	6	06-06-2024	Paragraaf 1, 1.8.4, 2.2, 2.4, 3.7, 4.6, 6.1 bijge- werkt, 4.6.6 toegevoegd.
2	7	05-12-2024	Paragraaf 1.8.4, 3.7, 4.6, 6.5, 7.4.1.2, 7.5.18, 10.1 bijgewerkt, 2.12 toegevoegd.

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

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



# Inhoudsopgave

1	Over	deze handleiding	8
	1.1	Toepassingsgebied van deze handleiding	8
	1.2	Fabrikant	8
	1.3	Beoogd gebruik	8
	1.4	Gebruiksbereik/toepassingsbereik	9
	1.5	Onjuist gebruik	9
	1.6	Garantie	9
	1.7	Handelsmerken	9
	1.8	Referentiedocumenten	9
	1.9	Conformiteit met wetten en normen	10
	1.10	Documentconventies	10
2	Veilig	heid	12
	2.1	Conventies voor veiligheidsmeldingen	12
	2.2	Algemene veiligheidsinformatie	14
	2.3	Privacyverklaring camera	16
	2.4	Toepassingsrisico's	16
	2.5	Exploitant	25
	2.6	Methode- en procesvalidatie	25
	2.7	Gebruikerskwalificatie	26
	2.8	Veiligheidselementen	27
	2.9	Productveiligheidstekens	33
	2.10	Laserstraling	38
	2.11	Ontsmettingsverklaring	39
	2.12	Incident melden	40
3	Techi	nische gegevens	41
	3.1	Typeplaatje	41
	3.2	Label met serienummer	42
	3.3	Afmetingen en gewichten	43



	3.4	Stroomvoorziening	44
	3.5	Gegevens- en stroomaansluitingen	45
	3.6	Omgevingsvoorwaarden	45
	3.7	Emissie en immuniteit	46
4	Besc	hrijving van de functie	48
	4.1	Overzicht	48
	4.2	Dek	48
	4.3	Robotarmen	52
	4.4	Vloeistofsysteem (Liquid FCA)	58
	4.5	Wassysteem (MCA 384)	59
	4.6	Opties en apparaten	59
_	Duri		
5	Bestl		71
	5.1	Bedieningselementen	/1
	5.2	Gebruikersinterface	72
	5.3	Foutsignalen en instrumentstatus	77
	5.4	Fluent ID-statusleds	79
6	Bedie	ening	80
Ŭ	6 1	Veiligheidsinstructies voor dit hoofdstuk	80
	6.2	Redriifem odi	81
	6.3	Inhadriifstalling	82
	6.4	Alverens con methodo to starton	02
	6.5	For methode uitvoron	90
	0.5		90
	0.0		107
	0.7		109
	0.8	Het Instrument uitschäkelen	
7	Syste	emverzorging	113
	7.1	Ontsmetting	113
	72	Beinigingsmiddelen	114
	1.2		
	7.3	Systeemverzorgingsmodus	115



	7.4	Systeemverzorgingstabellen	117
	7.5	Systeemverzorgingsactiviteiten	127
8	Verhe	elpen van storingen	157
	8.1	Veiligheidsinstructies voor dit hoofdstuk	157
	8.2	Tabellen voor het verhelpen van storingen	157
	8.3	Verhelpen van storingen handelingen	170
9	Verpa	akken, uitpakken, transport, opslag en verwijdering	200
9	Verpa 9.1	akken, uitpakken, transport, opslag en verwijdering Verpakkingslabels	<b>200</b> 200
9	Verpa 9.1 9.2	<b>akken, uitpakken, transport, opslag en verwijdering</b> Verpakkingslabels Afvoer	<b>200</b> 200 201
9	Verpa 9.1 9.2	<b>akken, uitpakken, transport, opslag en verwijdering</b> Verpakkingslabels Afvoer	<b>200</b> 200 201
9 10	Verpa 9.1 9.2 Klant	akken, uitpakken, transport, opslag en verwijdering Verpakkingslabels Afvoer	<ul><li>200</li><li>200</li><li>201</li><li>203</li></ul>
9 10	Verpa 9.1 9.2 Klant 10.1	akken, uitpakken, transport, opslag en verwijdering Verpakkingslabels Afvoer support Contacten	<ul><li>200</li><li>200</li><li>201</li><li>203</li><li>203</li></ul>
9 10	Verpa 9.1 9.2 Klant 10.1	akken, uitpakken, transport, opslag en verwijdering Verpakkingslabels Afvoer support Contacten	<ul><li>200</li><li>201</li><li>203</li></ul>

Inhoudsopgave





# 1 Over deze handleiding

Deze Bedieningshandleiding beschrijft de en alle informatie die nodig is om deze veilig te bedienen en in goede werkende staat te houden. Deze handleiding moet zorgvuldig worden gelezen, alvorens werkzaamheden uit te voeren aan de en deze te gebruiken.

Dit hoofdstuk beschrijft het doel van deze handleiding en specificeert het desbetreffende product. Verder wordt het gebruik van symbolen en conventies uitgelegd, evenals verdere algemene informatie.

Deze handleiding heeft betrekking op het -instrument zelf. Raadpleeg de specifieke handleidingen voor de modules voor belangrijke informatie over de submodules.



Deze Bedieningshandleiding bevat geen beschrijving van de software. Raadpleeg de bijbehorende softwarehandleiding voor meer informatie over de software. Zie paragraaf .

# 1.1 Toepassingsgebied van deze handleiding

Deze handleiding heeft betrekking op:

- Fluent 480 (onderdeelnummer 30042011)
- Fluent 780 (onderdeelnummer 30042021)
- Fluent 1080 (onderdeelnummer 30042031)

## **1.2 Fabrikant**

Adres van fabrikant



Tecan Schweiz AG Seestrasse 103 CH-8708 Männedorf Zwitserland

# 1.3 Beoogd gebruik

De Fluent is een volledig geautomatiseerd laboratoriumvloeistofverwerkingsplatform voor algemeen laboratoriumgebruik. Hij is bedoeld voor routinematige laboratoriumtaken, zoals pipetteren, vloeistofverwerking en robotgebaseerde manipulatie van laboratoriummateriaal bij bepaalde testprocedures.



### Voor dit beoogde gebruik is de Fluent Gx Assurance-software vereist.

Bepaalde opties van Tecan en apparaten van derden die met de Fluent kunnen worden gebruikt zijn uitsluitend voor onderzoekstaken bedoeld (research-use-only: RUO).

Als een RUO-optie of -apparaat in de Fluent is geïntegreerd, of als de Fluent Gx Assurance-software niet is geïnstalleerd, wijzigt het beoogde gebruik in:

De Fluent is een volledig geautomatiseerd laboratoriumvloeistofverwerkingsplatform voor onderzoek en industriële toepassingen. Hij is bedoeld voor routinematige laboratoriumtaken, zoals pipetteren, vloeistofverwerking en robotgebaseerde manipulatie van laboratoriummateriaal bij



bepaalde testprocedures (bijv. celtests, biochemische tests en verbindingenbeheer). Dit instrument is niet bedoeld voor gebruik in klinische diagnoseprocedures.

# 1.4 Gebruiksbereik/toepassingsbereik

De Fluent kan in vele verschillende laboratoriumomgevingen voor het beoogde doel worden gebruikt.

In elke omgeving is het individuele laboratorium verantwoordelijk voor de validatie van de Fluent in combinatie met de specifieke, te gebruiken vloeistoffen en het laboratoriummateriaal in de toepassingsworkflow of toepassingsmethode van het laboratorium.

# 1.5 Onjuist gebruik

Onjuist gebruik kan het veiligheidsconcept van de Fluent negatief beïnvloeden.

- De Fluent mag niet worden gebruikt met opties of componenten die niet zijn goedgekeurd door Tecan.
- De Fluent is niet explosieveilig en mag niet worden gemonteerd op plaatsen waar explosiegevaar bestaat.
- De Fluent mag niet worden gebruikt, als functionele veiligheidsvoorzieningen niet aanwezig zijn.

## 1.6 Garantie

De Fluent mag niet worden gebruikt met opties of componenten die niet zijn goedgekeurd door Tecan.

Het gebruik van niet-goedgekeurde componenten kan het veiligheidsconcept van de Fluent negatief beïnvloeden.

Door het gebruik van niet-goedgekeurde componenten vervalt elke veiligheidsgarantie en de conformiteit met nationale en internationale normen, zoals vereist voor NRTL-certificering, door EG-richtlijnen etc.

## 1.7 Handelsmerken

De productnamen, zowel geregistreerde als niet-geregistreerde handelsmerken, die in deze handleiding worden genoemd, worden uitsluitend gebruikt voor identificatiedoeleinden en blijven het exclusieve eigendom van hun respectievelijke eigenaars. Om redenen van eenvoud worden handelsmerksymbolen als <sup>®</sup> en <sup>™</sup> in deze handleiding niet herhaald.

# 1.8 Referentiedocumenten

Deze paragraaf bevat een lijst met documenten die nodig of praktisch kunnen zijn bij gebruik van de Fluent.

Onderstaande document-ID's zijn basisnummers. Daarom bevatten ze geen informatie over de taal, de documentversie of het medium (gegevensopslagmedium, papieren versie, downloadbaar bestand enz.) van het document.





Op basis van uw orderconfiguratie gelden eventueel ook de bedieningshandleidingen voor optionele uitrusting.

Controleer het toepassingsgebied van het bijbehorende document om er zeker van te zijn dat u de correcte versie heeft.

De document-ID verwijst niet naar bestelgegevens. Let bij het plaatsen van bestellingen op het nummer op de verpakking, cd-behuizing enz.

## 1.8.1 Instrumenthandleidingen

- Bedieningshandleiding Fluent<sup>®</sup> (doc-ID 399706)
- Referentiehandleiding Fluent® (doc-ID 399937)

## 1.8.2 Softwarehandleidingen

- Softwarehandleiding voor Tecan Sample Tracking Add-On (doc-ID 393933)
- Softwarehandleiding voor FluentControl-toepassing (doc-ID 399935)
- Softwarehandleiding voor Introspect-toepassing (doc-ID 400733)
- Softwarehandleiding voor MissionControl-toepassing (doc-ID 401940)

## 1.8.3 Handleidingen QC-kit

- Toepassingshandleiding QC-kit (doc-ID 397069)
- Handleiding voor toepassingssoftware QC-kit (doc-ID 397070)

## 1.8.4 Andere referentiedocumenten

- Bedieningshandleiding Fluent<sup>®</sup> Carousel (doc-ID 398350)
- HEPA-kap (doc-ID Caron 70072)
- Toepassingshandleiding Frida Reader™ (doc-ID 401882)
- Bedieningshandleiding Te-Shake™ (doc-ID 391496)
- Bedieningshandleiding Te-VacS<sup>™</sup> (doc-ID 391236)
- Bedieningshandleiding Fluent<sup>®</sup> Stacker (doc-ID 398658)
- Bedieningshandleiding MIO2 (doc-ID 394934)
- Bedieningshandleiding Resolvex i300 (doc-ID 402756)

## **1.9** Conformiteit met wetten en normen

De volgende verklaringen en certificeringen zijn van toepassing op Fluent:

- EG-conformiteitsverklaring met de geldende EU-richtlijnen (CE-markering)
- NRTL-certificering (National Recognized Testing Laboratory Nationaal erkend testlaboratorium)
- (IECEE) CB Scheme Certification (CB-markering)

Voor meer informatie over de markering, zie de paragraaf "Typeplaatje" [> 41].

## 1.10 Documentconventies

#### Verwijzingen

Verwijzingen verschijnen als volgt—bijv.:

Zie paragraaf "Veiligheid" [▶ 12]

• "Veiligheid" verwijst naar de bijbehorende paragraaftitel



	Het paginanummer staat tussen vierkante haakjes				
Voorwaarden	Voorwaarden verschijnen als volgt—bijv.:				
	✓ "Algemene veiligheidsinformatie" werd gelezen.				
Tips	Aanvullende tips verschijnen als volgt—bijv.:				
1	Voor veiligheidsconventies en -symbolen, zie hoofdstuk "Veiligheid" [▶ 12].				
Afbeeldingen	De afbeeldingen tonen eventueel componentversies die niet relevant zijn voor uw Fluent.				



# 2 Veiligheid

Dit hoofdstuk beschrijft het veiligheidsconcept van de Fluent, verschaft algemene regels voor correct gedrag en waarschuwingen met betrekking tot gevaren door gebruik van de Fluent.

## 2.1 Conventies voor veiligheidsmeldingen

## 2.1.1 Signaalwoorden

Tab. 1: Signaalwoorden

Signaalwoord	Betekenis
▲ GEVAAR	Duidt op een gevaarlijke situatie die, indien deze niet wordt voorkomen, leidt tot ernstig letsel of de dood.
	Duidt op een gevaarlijke situatie die, indien deze niet wordt voorkomen, kan leiden tot ernstig letsel of de dood.
	Duidt op een gevaarlijke situatie die, indien deze niet wordt voorkomen, kan leiden tot licht of ernstig letsel.
LET OP	Duidt op een niet-gevaarlijke situatie, die echter, indien deze niet wordt voorkomen, kan leiden tot schade of defecten aan de uitrusting, of onjuiste procesresultaten.



## 2.1.2 Veiligheidssymbolen





# 2.2 Algemene veiligheidsinformatie

## 

De Fluent is ontworpen en ingebouwd overeenkomstig de huidige stand van de techniek en de erkende technische veiligheidsregels. Toch kunnen er risico's ontstaan voor gebruikers, eigendommen en het milieu als de Fluent zonder de nodige zorg en aandacht wordt gebruikt.

De veiligheid van alle gebruikers en het personeel hangt af van de strikte inachtneming van deze veiligheidsinstructies en het bewustzijn met betrekking tot de veiligheidswaarschuwingen in deze handleiding.

- De volgende algemene veiligheidsinformatie in acht nemen.
- Deze handleiding moet altijd beschikbaar zijn voor alle personen die de erin beschreven taken uitvoeren.
- Gebruik altijd de voedingskabel die bij het instrument is geleverd.
- Gebruik de voedingskabel niet met andere producten.
- Wettelijke voorschriften zoals lokale, staats- en federale wetten met betrekking tot het gebruik of de toepassing van, de omgang met gevaarlijke materialen in verband met de Fluent moeten strikt worden opgevolgd.
- De exploitant is verantwoordelijk voor het opstellen van instructies in overeenstemming met de bedrijfsprocedures en de lokale wettelijke voorschriften. Deze instructies van de exploitant moeten strikt worden nageleefd.
- De correcte omgevingsvoorwaarden voor opslag en bedrijf in acht nemen.
- Structurele wijzigingen aan veiligheidsuitrustingen zijn verboden.
- Beschadigde veiligheidsuitrustingen moeten onmiddellijk worden vervangen zoals beschreven in deze handleiding.
- De Fluent mag op geen enkele wijziging worden gewijzigd zonder raadpleging en schriftelijke toestemming vooraf van Tecan. Geautoriseerde wijzigingen aan het systeem mogen alleen worden uitgevoerd door een technicus die is gecertificeerd voor reparatie en aanpassing van de Fluent. Tecan wijst elke claim als gevolg van ongeoorloofde wijzigingen af.
- Brandgevaar door onjuist gebruik van de Fluent. De Fluent mag niet worden gemonteerd op plaatsen waar explosiegevaar bestaat.
- · Brandgevaar door ontvlambare vloeistoffen of systeemvloeistof.
- Vermijd de vorming en ophoping van brandbare dampen.
- Chemische, biologische en radioactieve gevaren kunnen ontstaan door stoffen die worden gebruikt, monsters en reagentia die worden verwerkt door de Fluent (bijv. tijdens in- en uitladen). Hetzelfde geldt voor de verwijdering van afval.
  - Houd altijd rekening met mogelijke gevaren die van deze stoffen uitgaan.
  - Geschikte veiligheidskleding, veiligheidsbril, adembescherming en veiligheidshandschoenen gebruiken.
  - De omgang met deze stoffen en de verwijdering van afval kan onderworpen zijn aan de lokale, landelijke of federale wetgeving, of aan voorschriften met betrekking tot gezondheid, milieu of veiligheid. De desbetreffende bepalingen strikt in acht nemen.



- Elke verontreiniging moet onmiddellijk worden behandeld zoals in deze handleiding beschreven.
- De gebruiker is ervoor verantwoordelijke dat de Fluent altijd wordt bediend onder geschikte omstandigheden, en dat onderhouds-, service- en reparatiewerkzaamheden zorgvuldig, op tijd en alleen door geautoriseerd personeel worden uitgevoerd.
- Risico van niet-correcte meetresultaten. Nadat systeemverzorging of onderhoud is uitgevoerd, mag de bediening alleen worden voortgezet nadat is gecontroleerd of aan alleen voorwaarden voor correcte systeemwerking is voldaan.
- Gebruik uitsluitend aanbevolen verbruiksmiddelen binnen de vervaldatum en originele reserveonderdelen voor onderhoud aan en reparatie van het systeem om een correcte en betrouwbare werking te garanderen.
- Als de huid in contact komt met systeemvloeistof op het instrument, kan dit leiden tot letsel.
  - Draag altijd beschermende kleding overeenkomstig GLP.
- Zwaar voorwerp! Til het instrument niet op.
- Bedien het systeem niet zonder dektrays en deksegmenten.
- Dektrays vangen gemorste vloeistoffen op in het handmatige laadgebied van het dek. Het systeem moet worden gebruikt met zo veel mogelijk onder het dek gemonteerde dektrays, om alle gemorste vloeistoffen op te kunnen vangen. Gebruik het systeem niet zonder dektrays.
- Als geen residu-overdracht is toegestaan, wordt het gebruik van wegwerpnaalden met filters sterk aanbevolen.
- Mogelijke botsing. Plaats geen apparaten zonder Tecan-modelgegevens op het dek.
- Uitbreiding 300 is ontworpen voor een maximale belasting van 40 kg (88 lbs) en is alleen bedoeld voor gebruik met opties met een gewicht tot 40 kg (88 lbs).
- De Fluent wordt geleverd met een label met een veiligheidssymbool voor biologische gevaren dat door de gebruiker moet worden aangebracht bij gebruik van biologisch gevaarlijke stoffen. Breng het label aan op de voordeur op een plaats die zichtbaar is voor de gebruiker en handig is voor de toepassing. Zie paragraaf Productveiligheidstekens.
- De opties die op het werkblad van de worden gebruikt, kunnen sterke magnetische velden opwekken die de werking van medische hulpmiddelen geïmplanteerd bij of gedragen door een gebruiker, zoals pacemakers of insulinepompen, kunnen verstoren. De wordt geleverd met een veiligheidsaanduiding van sterke magnetische velden, die door de gebruiker op het deurtje aan de voorkant moet worden aangebracht op een plaats die duidelijk zichtbaar is voor de gebruiker en die geschikt is voor de toepassing in geval van gebruik van opties die sterke magnetische velden genereren.
- De ethernetkabel van de DeckCheck-camera's wordt geïnstalleerd door een technicus en moet te allen tijde op de Fluent-pc zijn geïnstalleerd (EMC). De Ethernet-interface mag niet worden verbonden met een netwerk.
- Alleen voor inwoners van Californië: Dit product kan u blootstellen aan chemische stoffen zoals lood. Deze stoffen zijn in de staat Californië bekend als veroorzakers van kanker en geboorteafwijkingen of andere schade aan de voortplanting. Voor meer informatie, zie www.P65Warnings.ca.gov/product.



## 2.3 Privacyverklaring camera

Het Fluent-systeem is voorzien van camera's op het frontprofiel aan de binnenzijde. De camera's zijn gericht op het voor- en achterdek. Zicht naar beneden door de zijwanden van acrylglas is mogelijk.

- De gebruiker is ervoor verantwoordelijk dat alle mensen in de ruimte erover geïnformeerd zijn dat er camera's actief zijn.
- De gebruiker is ervoor verantwoordelijk dat het personeel niet kan worden geïdentificeerd aan de hand van gemaakte foto's, bijvoorbeeld als het instrument naast (bij) een werkplek staat of als er achter- of zijpanelen zijn uitgesneden of als een acrylglazen paneel wordt gebruikt ter vervanging van de achterwand.

Systeem- functie/mo- dule	Mogelijke foutmodus	Potentieel ge- volg van de fout	Mogelijke/potenti- ële oorzaak	Etikettering of mitigatie
Systeem	Onvoldoen- de onder- houd	Veiligheid of gezondheid van gebruikers: po- tentiële veront- reiniging van instrument	Gebruiksfout: niet in acht nemen van de bedienings- handleiding of on- derhoudsinstruc- ties	De gebruiker dient de juiste ver- bruiksmiddelen te gebruiken en de instructies voor preventief on- derhoud te volgen (raadpleeg Systeemverzorging). De gebruiker dient beschermen- de kleding, handschoenen en een veiligheidsbril te dragen in overeenstemming met GLP en de geldende plaatselijke voorschrif- ten.
Systeem	Brand	Veiligheid of gezondheid van gebruikers: brand in ope- ratorlaboratori- um (instrument in brand)	Gas uit vluchtige ontvlambare vloeistoffen; rond- vliegende vonken van elektroplaat	Het instrument is niet explosie- veilig en de klant moet ervoor zorgen dat er geen hoge damp- concentratie aanwezig is (raad- pleeg Algemene veiligheidsinfor- matie).
Module FCA en Air FCA	Slijtage van het Z-as- mechanisme (bovenge- middelde gebruik)	Veiligheid of klinische toe- standen van monster: po- tentieel onjuiste Z-positionering in laboratori- ummateriaal	Bovengemiddeld gebruik van het apparaat in com- binatie met ge- bruik van weg- werpnaalden Hoog percentage doorprikstappen in toepassing	Het systeem informeert de ge- bruiker als de Z-assen 90% van de verwachte levensduur van de as hebben bereikt.

# 2.4 Toepassingsrisico's



Systeem- functie/mo- dule	Mogelijke foutmodus	Potentieel ge- volg van de fout	Mogelijke/potenti- ële oorzaak	Etikettering of mitigatie
Module FCA en Air FCA	Slijtage van het P-as- mechanisme (bovenge- middeld ge- bruik)	Veiligheid of klinische toe- standen van monster: po- tentieel onjuiste P-positionering in laboratori- ummateriaal	Bovengemiddeld gebruik van het apparaat in com- binatie met ge- bruik van weg- werpnaalden Hoog percentage doorprikstappen in toepassing	Het systeem informeert de ge- bruiker als de P-assen 90% van de verwachte levensduur van de as hebben bereikt.
Module FCA en Air FCA	Slijtage van X-aandrijf- tandwiel (bovenge- middeld ge- bruik)	Veiligheid of klinische toe- standen van monster: po- tentiële veront- reiniging van monsters met polyamidedeel- tjes	Bovengemiddeld gebruik van appa- raat in combinatie met het plaatsen van kritisch labo- ratoriummateriaal aan achterzijde van instrument	Vermijd het plaatsen van deel- tjesgevoelige elementen (bijv. monsters en reagentia) aan de achterzijde van het instrument of breng een deeltjesbescherming op de bovenzijde van het labora- toriummateriaal aan (bijv. dek- sels).
Module FCA en Air FCA	Storende signalen als gevolg van doorprikken van septum	Veiligheid of klinische om- standigheden van het pati- ëntmonster: onjuiste cLLD - dit leidt tot luchtopzuiging en mogelijk onjuiste resul- taten	Interactie van de naald met sep- tum/folie	Werk uitsluitend met niet-gelei- dende folies voor doorpriktoe- passingen in combinatie met vloeistofniveaudetectie op de FCA en Air FCA. Raadpleeg de referentiehandleiding. De gebruiker dient de vloeistof- detectie in combinatie met door- prikken voor FCA en Air FCA te valideren.
Module FCA en Air FCA	Onjuiste monsterbe- handeling, onjuiste cLLD door schuim of bellen in re- agensflacon	Procesveilig- heid: onjuist verwerkte mon- sters	Bellen of schuim in de reagensfla- con veroorzaken een onjuiste cLLD en mogelijke op- zuiging van lucht met FCA of Air FCA	De gebruiker is verantwoordelijk voor het valideren van de toe- passing/het proces met betrek- king tot de juiste monstervoorbe- reiding.
Module FCA en Air FCA	Naaldblok- kering	Veiligheid of klinische toe- standen van patiëntmonster: potentieel on- juist gepipet- teerd volume	Opzuiging aan onderzijde van well (naaldblokke- ring)	De gebruiker dient de toepassing te valideren en zo te voorkomen dat de opzuiging te veel het ni- veau Z-max van aangepast la- boratoriummateriaal nadert.



Systeem- functie/mo- dule	Mogelijke foutmodus	Potentieel ge- volg van de fout	Mogelijke/potenti- ële oorzaak	Etikettering of mitigatie
Module FCA en Air FCA	FCA-buis- systeem: groei van micro-orga- nismen	Veiligheid of klinische toe- stand van het patiëntmonster: onjuist gepi- petteerd volu- me of veront- reiniging van monsters	Groei van micro- organismen (bio- film op de bin- nenzijde)	Gebruik gedeïoniseerd water als systeemvloeistof voor FCA en voer dagelijks onderhoud uit om het systeem te spoelen volgens de instructies onder dagelijkse systeemverzorging (raadpleeg Systeemverzorging) en houd u daarbij aan de toegestane reini- gingsmiddelen voor die stap.
Module MCA 96 en/ of MCA 384/9 6	Overstromen van mon- stervloeistof in multititer- plaat tijdens pipetteren	Veiligheid of klinische toe- stand van pati- ëntmonster: Mogelijke kruisbesmetting van monsters (overstromen)	Onjuist gedefini- eerde Z-niveaus door gebruiker (bijv. opzuiging van positie Z- max)	Definieer veilige posities voor opzuiging en afgifte. Raadpleeg de referentiehandleiding.
Module MCA 96 en/ of MCA 384/9 6	Monsters missen ge- heel of ge- deeltelijk de beoogde positie in de modus vrij dispenseren	Veiligheid of klinische toe- stand van pati- ëntmonster: mogelijke kruisbesmetting	Elektrostatische ladingen aan het naalduiteinde door gebruik van het instrument buiten de gespecificeer- de voorwaarden leiden ertoe dat het monster aan de naald blijft hangen of dat er ongecontroleerd wordt gesproeid	De gebruiker dient zich te houden aan de gespecificeerde ge- bruiksvoorwaarden voor MCA- vloeistofverwerking, met name instructies over de minimaal ver- eiste luchtvochtigheid (zie "Om- gevingsvoorwaarden" [> 45]). De gebruiker dient de afgifte- hoogte altijd binnen de well in te stellen. Raadpleeg de referentie- handleiding.
Module MCA 96 en/ of MCA 384/9 6	Lucht men- gen in plaats van vloeistof (monster/ reagens) voor mix-pi- petteren	Veiligheid of klinische toe- stand van pati- ëntmonster: mogelijk on- juiste verwer- king van mon- sters leidt tot verkeerde re- sultaten	Niet-geschikte trackingparame- ters door onjuiste combinatie van naalden en multi- titerplaten	De gebruiker dient de echte en de virtuele werktafel te vergelijken met elkaar aan de hand van de naam van het laboratoriummate- riaal in de virtuele werktafel. De gebruiker dient op het unieke kleurontwerp (specifiek voor naaldtype) en etikettering (voor filter en niet-filter) van DiTi-do- zen te letten. De gebruiker dient de werktafel- indeling te controleren alvorens een proces te starten.



Systeem- functie/mo- dule	Mogelijke foutmodus	Potentieel ge- volg van de fout	Mogelijke/potenti- ële oorzaak	Etikettering of mitigatie
Specifiek voor weg- werpnaald	DiTi's opha- len: Onjuist naaldtype gemonteerd	Veiligheid of klinische toe- stand van pati- ëntmonster: mogelijk geen of te weinig monster opge- zogen mogelijk kruis- besmetting van monsters	Gebruiksfout: Onjuiste dekinde- ling: de gebruiker plaatst de naald- doos op de ver- keerde plaats: naalden zijn korter dan verwacht Onjuiste dekinde- ling: gebruiker plaatst naalddoos met ongefilterde naalden in plaats van gefilterde naalden op werk- tafel Onjuiste dekinde- ling: gebruiker plaatst naalddoos op verkeerde po- sitie: naald heeft kleiner volume dan verwacht (bijv. 100 µl in plaats van 200 µl); naaldlengte zoals verwacht; vloeistof opgezogen in MCH	De gebruiker dient de echte en de virtuele werktafel te vergelijken met elkaar aan de hand van de naam van het laboratoriummate- riaal in de virtuele werktafel. De gebruiker dient op het unieke kleurontwerp (specifiek voor naaldtype) en etikettering (voor filter en niet-filter) van DiTi-do- zen te letten. De gebruiker dient de werktafel- indeling te controleren alvorens een proces te starten. Het mechanische ontwerp ga- randeert de zichtbaarheid van het witte filter. De referentiehandleiding bevat informatie over kleurcodering van DiTi-dozen en het verschil in lengte, en gefilterde DiTi's. Raadpleeg de referentiehandlei- ding.
Specifiek voor weg- werpnaald	Niet alle naalden val- len: sommi- ge veront- reinigde naalden blij- ven aan de kop hangen en vallen op monsterpla- ten	Veiligheid of klinische toe- stand van pati- ëntmonster: mogelijke kruisbesmetting	Elektrostatische lading veroorzaakt	De gebruiker dient zich te houden aan de gespecificeerde ge- bruiksvoorwaarden voor MCA- vloeistofverwerking, met name instructies over de minimaal ver- eiste luchtvochtigheid (zie "Om- gevingsvoorwaarden" [> 45]). Wegwerpnaalden zijn niet be- doeld voor hergebruik.
Module RGA	Plaatverlies door botsen met niet- uitgelijnd la- boratorium- materiaal	Procesveilig- heid: plaatver- lies, verlies van monsters	Als meer dan 4 multititerplaten zijn gestapeld, kan dit tijdens transport leiden tot onjuiste uitlij- ning	Plaatbewegingen moeten worden gevalideerd voordat scripts met echte monsters worden uitge- voerd. Raadpleeg de controlelijst voor validatie in de software- handleiding van de toepassing.



Systeem- functie/mo- dule	Mogelijke foutmodus	Potentieel ge- volg van de fout	Mogelijke/potenti- ële oorzaak	Etikettering of mitigatie
Module Flu- entControl- software	Worktable- Base: on- juiste DiTi- status ge- rapporteerd	Procesveilig- heid: kruisbe- smetting/on- juiste resultaten	Kruisbesmetting door onjuiste in- formatie over ge- bruiksstatus van naalden	Gebruik 'Set Tips Back' (Naalden terugzetten) niet als foutmodus tot ernstige risico's leidt.
Module Flu- entControl- software	Core.Scrip- ting.Pro- gramming SetVariable bij runtime: onjuiste waarde	Procesveilig- heid: onjuiste resultaten	Fout in software: variabele is inge- steld op onjuiste waarde	Valideer de toepassing voor de specifieke bron, het doel en de bereiken van de variabele. Raad- pleeg de controlelijst voor valida- tie in de softwarehandleiding van de toepassing.
Module Flu- entControl- software	Core.Scrip- ting.Pro- gramming QueryVaria- ble bij runti- me of scriptstart: onjuiste UI- presentatie/ acceptatie van UI- waarde	Procesveilig- heid: onjuiste resultaten	Numerieke waarde wordt verkeerd geformatteerd of geconverteerd in UI	Valideer de toepassing voor de specifieke bron, het doel en de bereiken van de variabele. Raad- pleeg de controlelijst voor valida- tie in de softwarehandleiding van de toepassing.
Module Flu- entControl- software	Core.Scrip- ting.Pro- gramming ImportVaria- ble bij runti- me: onjuiste waarde ge- ïmporteerd	Procesveilig- heid: onjuiste resultaten	Onjuiste waarde wordt opgehaald van importbron	Valideer de toepassing voor de specifieke bron, het doel en de bereiken van de variabele. Raad- pleeg de controlelijst voor valida- tie in de softwarehandleiding van de toepassing.
Module Flu- entControl- software	Core.Scrip- ting.Pro- gramming ExportVaria- ble bij runti- me: onjuiste waarde naar bestand ge- ëxporteerd	Procesveilig- heid: onjuiste resultaten	Onjuiste waarde wordt naar ex- portbestand ge- schreven	Valideer de toepassing voor de specifieke bron, het doel en de bereiken van de variabele. Raad- pleeg de controlelijst voor valida- tie in de softwarehandleiding van de toepassing.



Systeem- functie/mo- dule	Mogelijke foutmodus	Potentieel ge- volg van de fout	Mogelijke/potenti- ële oorzaak	Etikettering of mitigatie
Module Flu- entControl- software	API: Varia- bele opha- len/instellen of expressie omzetten mislukt	Procesveilig- heid: onjuiste resultaten	Verkeerde varia- belewaarde opge- haald of verkeerde waarde toegewe- zen/verkeerd ex- pressieresultaat geretourneerd	Valideer de toepassing voor de specifieke bron, het doel en de bereiken van de variabele. Raad- pleeg de controlelijst voor valida- tie in de softwarehandleiding van de toepassing.
UVC-licht	Onjuist ge- bruik in toe- passing	Gebrek aan ef- fectiviteit	Onjuist gebruik in toepassing	Zie de specifieke instructies in paragraaf "Optische straling (UVC)" [> 32].
Buisrotator/ doorprik- naalden (Mix & Pierce)	Onjuist ge- bruik in toe- passing	Gebrek aan ef- fectiviteit	Onjuist gebruik in toepassing	Zie de specifieke instructies in paragraaf "Mix & Pierce" [> 63].
Frida Reader	Onjuist ge- bruik in toe- passing	Gebrek aan ef- fectiviteit	Onjuist gebruik in toepassing	Zie de specifieke instructies in paragraaf "Frida Reader" [> 66].
Alle	Ineffectief gebruik in toepassing	Gebrek aan ef- fectiviteit in toepassing	Gebrek aan sys- teemverzorging	Zie onderdeelspecifieke instruc- ties in het hoofdstuk Systeem- verzorging.
Verwerking van potenti- eel gevaar- lijke materi- alen	Verontreini- ging met potentieel gevaarlijke materialen	Mogelijke ge- varen voor ge- bruikers, ei- gendom en het milieu.	Niet-naleving van algemene veilig- heidsinformatie	Raadpleeg de onderdeelspecifie- ke instructies in paragraaf Alge- mene veiligheidsinformatie.
MCA 96, cLLD	Onjuiste meetresul- taten	Onjuiste me- ting: het gede- tecteerde vloei- stofniveau geldt niet voor alle wells van de multititer- plaat: onjuist testresultaat of verlies van het monster	Gebruik van onge- schikt/incompati- bel laboratorium- materiaal	De cLLD-functie kan dan alleen worden gebruikt in trays.



Systeem- functie/mo- dule	Mogelijke foutmodus	Potentieel ge- volg van de fout	Mogelijke/potenti- ële oorzaak	Etikettering of mitigatie
MCA 96, al- gemeen	Incompati- biliteit van verbruiks- middelen/ onderdelen/ modules	DiTi's verloren tijdens proces: mogelijk on- herstelbaar verlies van het monster. Mo- gelijke kruisbe- smetting.	Gebruik van onge- schikte/incompa- tibele verbruiks- middelen. Er wordt een op- dracht gegeven om een offset op te pakken die niet compatibel is met de doos of tray waaruit de DiTi's moeten worden geplaatst (bijv. onjuiste versie van DiTi-tray). DiTi's zijn onjuist ge- plaatst en zijn ver- loren gegaan tij- dens het proces	Gebruik DiTi-dozen met Tecan- combitray.
MCA 96, cLLD	Onjuiste meetresul- taten	Onjuiste me- ting: Fout-po- sitieve detectie van cLLD-sub- systeem.	De gebruiker plaatst andere Di- Ti's op de werkta- fel dan in de soft- ware wordt weer- gegeven. Het aantal geplaatste DiTi's wijkt af van het aantal DiTi's dat door de soft- ware wordt ver- wacht als gevolg van een handeling van de gebruiker (de gebruiker ver- wijdert bijvoor- beeld enkele Di- Ti's). Fout-posi- tieve detecties als gevolg van onjuis- te cLLD-drempel.	Belangrijk: het aantal DiTi's dat wordt gebruikt voor cLLD-detec- tie, moet overeenkomen met het aantal dat wordt weergegeven in de software.



Systeem- functie/mo- dule	Mogelijke foutmodus	Potentieel ge- volg van de fout	Mogelijke/potenti- ële oorzaak	Etikettering of mitigatie
MCA 96, al- gemeen	Verontreini- ging van monster	Verontreiniging van monster als gevolg van morsen na een botsing. Onjuist testresultaat of verlies van het monster.	Onjuiste parame- ters gebruikt (on- juiste vector voor oppakken of neerzetten, onjuist type gereed- schap). Botsing van de kop van de MCA 96 tijdens het verplaatsen van de vector. Bij- voorbeeld met hoog laboratori- ummateriaal op aangrenzende nesten.	Er zijn risico's verbonden aan onjuist gebruik van de vectorver- plaatsingsfunctie.
MCA 96, al- gemeen	Verontreini- ging van monster	Veiligheid of klinische toe- stand van pati- ëntmonster: Verontreiniging van monster: mogelijke kruisbesmetting door slijtage van de grijper- band.	Slijtage van band- materiaal gedu- rende de levens- duur: Door slijtage kunnen stukjes van de band van de G-as in pati- ëntmonsters op de werktafel te- rechtkomen en uiteindelijk het monster chemisch verontreinigen.	Door slijtage kunnen stukjes (van banden) en stof op de werktafel terechtkomen en monsters/che- micaliën vervuilen. Om dit te voorkomen, moet laboratorium- materiaal dat dergelijke gevoelige vloeistoffen bevat, worden be- schermd door deksels te plaat- sen.
MCA 96, al- gemeen	Verontreini- ging van monster	Veiligheid of klinische toe- stand van pati- ëntmonster: Mogelijk on- juiste resulta- ten: mogelijke kruisbesmetting door morsen in aangrenzende plaatholten.	Spatten/morsen van monstervloei- stof. Verontreini- ging van DiTi-ke- gels door vloeistof in vastgehouden laboratoriumma- teriaal, bijvoor- beeld wanneer volledig laborato- riummateriaal bij het afval wordt gedeponeerd en er vloeistof wordt gemorst.	Maak het laboratoriummateriaal eerst leeg voordat u dit afvoert.



Systeem- functie/mo- dule	Mogelijke foutmodus	Potentieel ge- volg van de fout	Mogelijke/potenti- ële oorzaak	Etikettering of mitigatie
MCA 96, al- gemeen	Residu- overdracht van monster of reagens (FC)	Verontreinigde DiTi-kegels en cilinderblokken, wat kan leiden tot mogelijk kruisbesmette monsters.	Het apparaat pro- beert DiTi's op te pakken. Aange- zien de software geen ontbrekende DiTi's kan detec- teren, gaat het apparaat verder zonder foutmel- ding. De DiTi-ke- gels komen mo- gelijk in contact met de bak en zuigen vloeistof op uit de bak.	De MCA 96 kan geen lege naalddozen detecteren bij ge- bruik van DiTi-doostypen waarbij de rand van de naald gelijk ligt met de bovenkant van de doos.
MCA 96, al- gemeen	Spatten/ morsen van monster- vloeistof	Verontreiniging van het mon- ster als gevolg van morsen vanwege on- voldoende ste- vig laboratori- ummateriaal.	Spatten/morsen van monstervloei- stof door gebruik van niet stevig ge- noeg laboratori- ummateriaal.	Om vervorming van laboratori- ummateriaal te voorkomen, mag u uitsluitend stevig genoeg labo- ratoriummateriaal gebruiken om vloeistoffen te transporteren.
MCA 96, al- gemeen	Grijpervinger kan bescha- digd raken na een bot- sing	Grijpervinger beschadigd		Controleer de grijpervingers en vervang ze indien beschadigd.
Mix & Pier- ce-wassta- tion	Verontreini- ging van monster	Het afval werd niet goed af- gevoerd. Verontreiniging van het mon- ster door on- juist drainen van het afval en onjuist was- sen van de naalden.	Onjuist gewassen naalden door overstroming van het wasstation veroorzaakt door verstopping van de aansluiting op het wasstation.	Het systeem moet regelmatig worden onderhouden. De connectoren van het wassta- tion moeten binnen cycli van 2- 3 maanden worden vervangen voor volbloedtoepassingen om verstopping te voorkomen. Om corrosie te voorkomen, mo- gen potentieel corrosieve vloei- stoffen, zoals 2% bleekmiddel, niet worden afgevoerd via het wasstation, de connectors en de slangen zonder extra te spoelen met neutrale vloeistoffen, zoals water.



# 2.5 Exploitant

De exploitant moet ervoor zorgen dat de Fluent, en in het bijzonder de veiligheidskenmerken, goed functioneren en dat al het personeel dat met het instrument in contact komt voldoende is geïnstrueerd.

## Verantwoordelij

kheden

- Methode- en procesvalidatie.
- Definiëren van de processen in overeenstemming met de standaard werkprocedures.
- Ervoor zorgen dat de installatie- en operationele kwalificaties (IQ OQ's) zijn afgesloten.
- Zorg ervoor dat al het personeel dat met de Fluent in contact komt adequaat is geïnstrueerd.
- Ervoor zorgen dat geschikte beschermende kleding en beschermingsmiddelen beschikbaar zijn.
- Ervoor zorgen dat het onderhoud is uitgevoerd en de Fluent veilig wordt gebruikt.
- Ervoor zorgen dat aan alle veiligheidsvoorschriften en -richtlijnen voor laboratoria wordt voldaan.

## 2.6 Methode- en procesvalidatie

Neem bij het valideren van de methode en het proces het volgende in acht:

- Bij gebruik van vaste naalden met MCA 384 of FCA moet u ervoor zorgen dat de wasprocedure effectief is voor het verwachte concentratiebereik van het monster en de gevoeligheid van de test.
- Controleer of de gepipetteerde volumes voldoen aan de precisie- en nauwkeurigheidseisen van het geautomatiseerde proces.
- Bij gebruik van niet-Tecan of aangepast laboratoriummateriaal en bij het opzuigen met tracking moet ervoor worden gezorgd dat de reservoirdefinitie correct is (bijv. dat de juiste snelheid voor tracking wordt gebruikt) om het opzuigen van lucht te vermijden.
- De functie van de fasescheider is geverifieerd voor gebruik met standaard
  Tecan-wegwerpnaalden van 1 ml en Tecan-wegwerpnaalden met brede
  opening van 1 ml. Raadpleeg de referentiehandleiding (zie
  "Referentiedocumenten" [> 9]) voor meer informatie over ondersteunde Tecan-verbruiksmiddelen.
- Verantwoordelij kheid van de key-operator
- Valideer de vloeistofdetectie op het Fluent Stacker-transferstation.
- Valideer het correcte gebruik van het MCA-wasstation door de toepassing.
- Valideer de toepassing met betrekking tot correcte pipetteervolumes en tracking.
- Valideer de toepassing om opzuiging te dicht bij Z-max van aangepast laboratoriummateriaal te voorkomen.
- Valideer doorpriktoepassingen met betrekking tot de vereiste neerhouders (actief of passief).
- Als chemicaliën en laboratoriummateriaal niet worden verwijderd, moet het effect van UVC-licht op chemicaliën en laboratoriummateriaal op het dek worden geëvalueerd en moet de test worden gevalideerd.
- Voer na de cyclus een handmatige controle op correcte pipetteervolumes uit.



 Personeel moet worden geïnformeerd over de privacyverklaring bij cameragebruik (zie "Privacyverklaring camera" [> 16]).

# 2.7 Gebruikerskwalificatie

Het laboratoriumpersoneel moet volledig gekwalificeerd en opgeleid zijn voor het werken met de Fluent. De werkzaamheden die in dit Bedieningshandleiding worden beschreven mogen alleen worden uitgevoerd door bevoegd personeel met de hieronder voorgeschreven kwalificaties.

Laboratoriumpersoneel moet:

- een geschikte technische opleiding hebben,
- vertrouwd zijn met de laboratorium-veiligheidsregels en richtlijnen,
- vertrouwd zijn met de instructies voor de veiligheidselementen van het instrument,
- beschermende kleding en beschermingsmiddelen gebruiken,
- vertrouwd zijn met en zich houden aan goede laboratoriumpraktijken,
- en de instructies in de handleiding hebben gelezen en begrepen.

Tecan adviseert dat de operator deelneemt aan een operatoropleiding. Raadpleeg de Tecan-klantenservice over beschikbare opleidingen. Zie paragraaf "Klantsupport" [▶ 203].

#### 2.7.1 Operator

De operator (laboratoriumtechnicus) werkt voor de exploitant.

- Geen specifieke toepassing of systeemkennis
  - Beheersing van de ter plaatse gebruikte talen
  - Beheersing van Engels wordt aanbevolen

De operator heeft toegangsrechten voor de toepassingssoftware waarmee hij methoden en systeemonderhoud kan uitvoeren en de benodigde training van de key-operator ontvangt.

#### 2.7.2 Key-operator

•

De key-operator (toepassingsspecialist) ondersteunt de exploitant of werkt voor dezelfde onderneming.

Vereiste vaardigheden Verantwoordelij

Vereiste

vaardigheden

- Uitgebreide kennis van de toepassingBegrensde systeemkennis
  - Beheersing van de ter plaatse gebruikte talen
  - Beheersing van het Engels

Instrueren van de operator

• Diepgaande kennis van de bijbehorende softwarehandleiding

Verantwoordelij kheden

- Schrijven, uitvoeren en valideren van methodes
  - Helpen van de operator bij het oplossen van problemen met het instrument



# 2.8 Veiligheidselementen

## **⚠ VOORZICHTIG**

#### Bewegende delen

De beschermings- en veiligheidsvoorzieningen op de Fluent mogen tijdens het gebruik niet verwijderd, gedeactiveerd of overbrugd worden.

 Als er apparaten worden verwijderd (bijv. voor onderhoudswerkzaamheden), moeten alle beschermings- en veiligheidsvoorzieningen opnieuw geïnstalleerd, geactiveerd en gecontroleerd worden voordat het gebruik kan worden voortgezet.

Veiligheidspanelen en veiligheidssensors zijn integrale delen van de Fluent, terwijl vergrendelingen van deuren en kabinetten alleen bij bepaalde systemen aanwezig kunnen zijn.

## 2.8.1 Veiligheidspanelen

Fluent is beschermd met veiligheidspanelen:

De **veiligheidspaneel aan de voorzijde** kan worden geopend en is uitgerust met deursensors die een actieve stop activeren. Het veiligheidspaneel aan de voorzijde kan worden afgesloten met optionele deurvergrendelingen.

De Fluent in combinatie met MCA 96- of MCA 384-armen mag alleen worden gebruikt met het volledige veiligheidspaneel aan de voorzijde.

De klep wordt ondersteund door gasveren. De operator moet de klep volledig openen voor veilige/volledige toegang tot het instrument.

De **verdunnerpaneel** kan worden geopend zonder de werking van Fluent te beïnvloeden (behalve voor Fluents met geïnstalleerde UVC-lichtoptie, de sensor van het verdunnerpaneel activeert een snelle stop wanneer het paneel wordt geopend).

De boven- en zijpanelen zijn vast bevestigd.

#### 2.8.1.1 Veiligheidspanelen aan de voorzijde

Het veiligheidspaneel aan de voorzijde voorkomt directe toegang tot de robotarmen en tot de elementen op het instrumentdek tijdens gebruik. Dit dient de veiligheid van het personeel en verbetert de veiligheid van de methode. Bovendien beschermt het veiligheidspaneel aan de voorzijde de gebruiker tegen het morsen van monsters of reagens. Er bestaan verschillende soorten veiligheidspanelen aan de voorzijde. **2 - Veiligheid** Veiligheidselementen



Volledig veiligheidspane el aan de voorzijde



Afb. 1: Volledig veiligheidspaneel aan de voorzijde

Het volledige veiligheidspaneel aan de voorzijde heeft de volgende kenmerken:

- Geen bewegende delen (bewegende delen, mechanische gevaren)
- · Bescherming van de monsters tegen externe invloeden (methodeveiligheid)
- Bescherming tegen morsen van monster of reagens



*Met volledige veiligheidspanelen aan de voorzijde is laden alleen per batch mogelijk.* 



Volledig

el aan de voorzijde (UVC)

veiligheidspane

Afb. 2: Volledig veiligheidspaneel aan de voorzijde (UVC)

Het volledige veiligheidspaneel aan de voorzijde (UVC) heeft de volgende kenmerken:

- Geen bewegende delen (bewegende delen, mechanische gevaren)
- Bescherming van de monsters tegen externe invloeden (methodeveiligheid)
- · Bescherming tegen morsen van monster of reagens
- Bescherming tegen optische straling (UVC)



Met volledige veiligheidspanelen aan de voorzijde is laden alleen per batch mogelijk.

# **A VOORZICHTIG**

#### Bewegende delen!

Bewegende MCA, FCA en Air FCA kunnen leiden tot handletsel, als er tijdens een cyclus door het halve veiligheidspaneel aan de voorzijde of door het veiligheidspaneel aan de voorzijde met uitbreiding in het instrument wordt gegrepen.

Grijp nooit tijdens een cyclus in het instrument.

2 - Veiligheid Veiligheidselementen





Half veiligheidspane el aan de voorzijde

Afb. 3: Half veiligheidspaneel aan de voorzijde

Het halve veiligheidspaneel aan de voorzijde heeft de volgende kenmerken:

- Beperkte toegang tot bewegende delen (bewegende delen, mechanische gevaren)
- · Bescherming tegen morsen van monster of reagens



Door het halve veiligheidspaneel aan de voorzijde heeft de operator slechts beperkte toegang tot het instrumentdek. Het in- en uitladen van runners is mogelijk zonder het paneel te openen; de operator kan monsters of reagentia herladen tijdens een methodecyclus.



Veiligheidspane el aan de voorzijde met uitbreiding

Afb. 4: Veiligheidspaneel aan de voorzijde met uitbreiding

Het veiligheidspaneel aan de voorzijde heeft de volgende kenmerken:

- Beperkte toegang tot bewegende delen (bewegende delen, mechanische gevaren)
- Bescherming tegen morsen van monster of reagens
- Maakt het gebruik mogelijk van een DiTi-afvoerstation aan de voorzijde dat uit het dek steekt en een omlaag gerichte opening in het veiligheidspaneel aan de voorzijde vereist.



Met het veiligheidspaneel aan de voorzijde met uitbreiding is laden alleen per batch mogelijk.

#### 2.8.1.2 Veiligheidspanelen voor optionele apparaten

Als een optioneel apparaat is toegevoegd aan of verwijderd van de zijkant van de Fluent, moet een geschikt zijveiligheidspaneel geïnstalleerd zijn. Raadpleeg de "Klantsupport" [> 203].

### 2.8.2 Instrumentdeurvergrendelingen (optioneel)

Twee optionele deurvergrendelingen kunnen voorkomen dat het veiligheidspaneel aan de voorzijde opengaat en beveiligen zo het actieve proces. Dit voorkomt een ongeoorloofde onderbreking van de procescyclus. Om het proces te stoppen kan op het touchscreen een pauze worden aangevraagd.



#### 2.8.3 Kabinetdeurvergrendelingen

Als een RGA-lange as onder het dek toegang heeft, moet de kabinetdeur die zich het dichtst bij het toegangspunt bevindt voorzien zijn van een optie voor een deurvergrendelingssensor. Als er meer dan één toegangspunt onder het dek is geïmplementeerd of als het toegangspunt tijdens de levensduur van het instrument werd gewijzigd, dan moet elke deur in de buurt van het toegangspunt worden uitgerust met een deurvergrendelingssensor.

Als het instrument een HEPA-kap heeft, moeten alle kabinetdeuren zijn uitgerust met een deurvergrendelingssensor.

#### 2.8.4 Optische straling (UVC)

De Fluent kan worden uitgerust met een optionele HEPA-kap met een UVC-licht of een afzonderlijke UVC-lichtoptie.

Blootstelling aan UVC-licht moet worden voorkomen omdat het kan leiden tot letsel. Het UVC-licht wordt automatisch uitgeschakeld als het veiligheidspaneel aan de voorzijde wordt geopend, en, bij de UVC-lichtoptie, ook als de verdunnerafdekking wordt geopend. Speciale UVC-resistente veiligheidspanelen worden op de Fluent met UVC-licht geïnstalleerd.

UVC-licht kan in ontsmettingsprocedures worden gebruikt. De geschiktheid en effectiviteit van UVC voor afzonderlijke processen moet door de gebruiker worden gevalideerd.



Zie ook de handleiding van de fabrikant van de HEPA-kap.

## 2.8.5 Externe deurvergrendelingen

Externe deurvergrendelingen worden aangebracht op Fluent-installaties in een externe behuizing. De deurpanelen van de externe behuizing vervangen de mechanische veiligheidsfunctie van het veiligheidspaneel aan de voorzijde en de kastdeuren van de Fluent en de externe deurdocks met geïntegreerde sensoren vervangen de deursensor- en deurvergrendelingsfuncties van het veiligheidspaneel aan de voorzijde en de kastdeuren van de Fluent.



Externe deurvergrendelingen laten geen ActiveStop toe. Een stop of pauze van het proces kan op het touchscreen worden aangevraagd.



# 2.9 Productveiligheidstekens

Om veiligheidsredenen zijn er op de Fluent veiligheidstekens aangebracht. Beschadigde, verloren of onleesbare veiligheidstekens moeten onmiddellijk zoals afgebeeld worden vervangen. Voor de betekenis van de veiligheidssymbolen, zie de paragraaf "Conventies voor veiligheidsmeldingen" [> 12].





Afb. 5: Standaardinstrument



Afb. 6: Instrument met UVC

UVC





Afb. 8: Biologisch gevaar





De Fluent wordt geleverd met een label met een veiligheidssymbool voor biologische gevaren dat door de gebruiker moet worden aangebracht bij gebruik van biologisch gevaarlijke stoffen.

Breng het label aan op de voordeur op een plaats die zichtbaar is voor de gebruiker en handig is voor de toepassing.



Afb. 9: Instrument met half veiligheidspaneel aan de voorzijde

Instrument met half veiligheidspane el aan de voorzijde Instrument met veiligheidspane el aan de

voorzijde met uitbreiding





Afb. 10: Instrument met veiligheidspaneel aan de voorzijde met uitbreiding

MCA



Afb. 11: Veiligheidssymbool op MCA 384


Dekuitbreiding



Afb. 12: Dekuitbreiding

#### 2.9.1 Mix & Pierce-werkstation



FCA beschermafdek king

Afb. 13: Beschermafdekking



# 2.10 Laserstraling

Fluent kan worden uitgerust met laserbarcodescanners. De laserstraling van deze barcodescanners is een gecollimeerde bundel in het zichtbare spectrum met een laag vermogen. De laserklassen van elke barcodescanner en van het gehele Fluent-systeem staan vermeld op het laserveiligheidslabel dat is aangebracht op de bijbehorende hardware.

Alle modules met lasers zijn geïdentificeerd met de desbetreffende laserveiligheidslabels.

Het Fluent instrument werd getest en gecertificeerd conform IEC 60825-1:2007 en IEC 60825-1:2014.



# 

De Fluent een klasse 1 laserproduct conform IEC 60825-1:2014 dat laserstraling uitzendt.

Verblinding, flitsblindheid en nabeelden kunnen door de laserstraal worden veroorzaakt.

• Staar niet in de laserstraal of in de reflecties ervan.

#### 2.10.1 Laserstralingsapparaten

Een standalone-barcodescanner kan op een apparaat worden gemonteerd.

Zorg ervoor dat het veiligheidslabel te allen tijde correct op de barcodescanner is aangebracht:

- Toelichting laserstralingslabel (A): Identificeert een KLASSE 2 LASERPRODUCT conform IEC 60825-1 met een geïntegreerd zichtbare, laserbarcodescanner met laag vermogen. Instrueert de gebruiker om niet in de laserstraal of de reflectie ervan te kijken.
- Klasse 2 lasers worden alleen gebruikt, als het systeem actief is en geen direct contact met de operator hebben.

Labelpositionering	Toelichting
A LASER RADIATION DO NOT STARE INTO BEAM Dutput 85,00 Multe duration 112,02 CLASS 2 LASER PRODUCT IEC/EN 60825-1 : 2007	Standalone-barcodescanner gemon- teerd op een hotel: label gepositioneerd onder de scan- ner.
LASER RADIATION Do Not & TARE INTO BEAM Output & 650m Wavelength & 650m Used to 122gs CLASS 2 LASER PRODUCT EC/EN 60825-1: 2007	Standalone-barcodescanner gemon- teerd op robotgrijperarm: label gepositioneerd op de scanner.



Labelpositionering	Toelichting
	Lading-ID: label gepositioneerd aan de achter- zijde van de scannerbehuizing.
AND EXPOSINE - LASE RADATOR	Lading-ID: label gepositioneerd aan de zijkant van de scannerbehuizing.

# 2.11 Ontsmettingsverklaring

Naast de regelmatige systeemverzorging en overeenkomstig de standaard laboratoriumvoorschriften moeten de Fluent en zijn onderdelen en accessoires onder de volgende omstandigheden grondig worden ontsmet:

- Voordat er aan de Fluent onderhouds- of servicewerkzaamheden worden uitgevoerd, in het bijzonder bij werkzaamheden door een FSE aan de Fluent
- Bij ongevallen (bijv. botsing, morsen etc.)
- Voordat de Fluent of delen of accessoires ervan naar Tecan (bijv. ter reparatie) worden teruggestuurd
- Voor opslag
- Voor uiteindelijke verwijdering
- In het algemeen voordat de Fluent of delen ervan worden verplaatst

De eigenaar van het instrument is volledig verantwoordelijk voor de effectieve ontsmetting van alle apparatuur.

Voor elke ingreep aan de Fluent door een technicus en voor het retourneren van de Fluent of onderdelen of accessoires ervan naar Tecan moet de eigenaar van het instrument het formulier voor de ontsmettingsverklaring invullen en ondertekenen om te bevestigen dat de ontsmetting is uitgevoerd in overeenstemming met de richtlijnen voor goede laboratoriumpraktijken. Neem contact op met uw lokale organisatie om dit formulier aan te vragen en lees paragraaf Ontsmetting.



Tecan behoudt zich het recht om de Fluent of onderdelen of accessoires ervan te weigeren of deze te bewerken, als er geen ontsmettingsverklaring wordt bijgevoegd.



# 2.12 Incident melden

Elk ernstig incident dat zich in verband met het hulpmiddel heeft voorgedaan, moet worden gemeld aan de fabrikant en de bevoegde autoriteit van de lidstaat waarin de gebruiker en/of de patiënt gevestigd is. Raadpleeg de paragraaf Fabrikant voor het adres van de fabrikant.



#### **Technische gegevens** 3

# 3.1 Typeplaatje



- Vor Service oder Wartungsarbeiten Netzstecker zienen
  Prior to any repair or maintenance job disconnect mains power

  - 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



When Laser Module(s) included VICLAS 11. 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"

 $(\epsilon)$ MET CAN ICES-001(B) / NMB-001(B)



Afb. 14: Typeplaatje

Het typeplaatje bevindt zich aan de achterzijde van de Fluent en bevat de volgende informatie:

Identificatiegegevens	Model	
	REF: Bestelgegevens (materiaalnummer)/revisieni- veau	
	Productiedatum (YYYYMMDD)	
	SN: Serienummer	
Technische gegevens	U, f: Voedingsspanning (volt), frequentie (Hertz)	
	P: Stroomverbruik (W)	
Adresgegevens	Naam en adres van de fabrikant	
Conformiteitsgege- vens	Conformiteitsmarkering	



# 3.2 Label met serienummer

Model	Instrument	Fluent 780	)	
REF	30042021	00		
	ECAN S H-8708	chwei Männe	iz AG edorf	
30	042021 00991	2123456		
<sup>SN</sup> 99	9121	234	56	

Afb. 15: Label met serienummer

Een serienummerlabel bevindt zich in de behuizing rechtsachter op het instrument en bevat de volgende gegevens:

Identificatiegegevens	Model
	REF: Bestelgegevens (materiaalnummer)/revisieni- veau
	SN: Serienummer
Adresgegevens	Naam en adres van de fabrikant





# 3.3 Afmetingen en gewichten

	Afmeting	Fluent 480	Fluent 780	Fluent 1080
а	Totale lengte	1150 mm (45,28 in.)	1650 mm (64,96 in.)	2150 mm (84,65 in.)
b	Voetafdrukdiepte	7	'80 mm (30,71 in	.)
С	Totale diepte	9	023 mm (36,34 in	.)
d	Totale hoogte op kabinet	1	977 mm (77,8 in	.)

Component	Fluent 480	Fluent 780	Fluent 1080
Basiseenheid	120 kg (264,5 lb.)	140 kg (308,6 lb.)	190 kg (418,9 lb.)
Verpakking	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.)		
384-kanaalkop	7,2 kg (15,9 lb.)		
RGA	10,2 kg (22,4 lb.)		
RGA-Z	10,6 kg (23,4 lb.)		



Component	Fluent 480	Fluent 780	Fluent 1080
сХР		1,2 kg (2,6 lb.)	
MCA 96 inclusief optionele grijper		19,7 kg (43,43 lb.)	

# 3.4 Stroomvoorziening

# LET OP

#### Oververhitting van de stroomvoorziening

De stroomvoorzieningseenheid kan beschadigd of onbruikbaar zijn.

- De stroomvoorziening mag niet worden afgedekt.
- De warmte van de stroomvoorziening moet kunnen worden afgevoerd.



Externe apparaten mogen niet op de stroomvoorziening worden aangesloten. Ze kunnen leiden tot een reset of stilstand van de Fluent

Tab. 2: Fluent stroomingang

Toevoer	Nominaal
Leidingspanning (enkele fase)	100-240 VAC
Ingangsstroom	9,8 A (bij 100 V) – 4 A (bij 240 V)
Frequentie	50–60 Hz

#### Tab. 3: Fluent stroomuitgang

Toevoer	Nominaal
Uitgangsspanning	24-28 V fabrieksinstelling: 25,2 V
Continustroom	500 W
Piekvermogen (begrensde tijd)	1500 W gedurende 3 seconden
Gewicht	3,8 kg (8,5 lbs.)

Max. spanningsschommeling van de netvoeding: ±10% van nominale spanning. Classificatie met betrekking tot de elektrische veiligheid volgens EN/IEC-normen:

#### *Tab. 4:* Elektrische specificaties (veiligheid)

Overspanning categorie	II	IEC 60664-1
Verontreinigingsgraad	2	(EN) IEC 61010-1





# 3.5 Gegevens- en stroomaansluitingen

Afb. 16: Gegevens- en stroomaansluitingen

- A Fluent instrument
- C Stroomvoorzieningseenheid
- I1 USB-interface

- Besturings-pc
- Wanduitgang
- Stroomkabel

De afbeelding toont de componenten van een monstersysteem met gegevens- en stroomaansluitingen. De Fluent-instrumentonderdelen zijn afgebeeld binnen de rechthoek. De hoofdschakelaar van het instrument maakt deel uit van de stroomvoorzieningseenheid. De stroomkabel is verbonden met een wanduitgang voor de netstroomvoorziening.

В

D

12

Gegevensoverdracht naar en van de Fluent wordt uitgevoerd via de USB-interface. De USB-kabel is verbonden met de pc die het instrument bestuurt.

# 3.6 Omgevingsvoorwaarden

# **⚠ VOORZICHTIG**

#### **Onjuiste pipetteervolumes**

Pipetteerresultaten kunnen door gebruiksomstandigheden worden beïnvloed.

Condensatie kan elektronische componenten beïnvloeden.

Als de Fluent wordt opgeslagen of getransporteerd bij temperaturen onder kamertemperatuur, moet hij na installatie enkele uren acclimatiseren.





S

Bedrijfsconditie

#### De Fluent mag uitsluitend binnen worden gebruikt en opgeslagen.

Bedrijfstemperatuur	15-32°C (59-90°F)
Bedrijfsvochtigheid	30-80% relatief (niet-condenserend) bij 30 °C (86 °F)
Bedrijfshoogte	max. 2000 m boven de zeespiegel

Bedrijfscondities voor vloeistofverwerking en pipetteren:

Ruimtetemperatuur	20-25°C (68-77°F)
Bedrijfsvochtigheid	30-60% relatief (niet-condenserend)
Bedrijfshoogte	max. 500 m boven de zeespiegel
Verdamping	In omgevingen met een verhoogde luchtstroom (door laminaire stro- ming, airconditioning of ventilatie etc.) wordt het risico op verdamping ook hoger. Dit kan de precisie van het pipetteren verminderen, in het bijzonder bij lage volumes of vluchti- ge stoffen.
	LET OP! Zorg ervoor dat de validatie condities overeenkomen met de cy- cluscondities.
Transporttemperatuur	-20 tot 60 °C (-4 tot 140 °F)
Transportvochtigheid	20-80% relatief (niet-condenserend
Opslagtemperatuur	1-60°C (34-140°F)
Opslagvochtigheid	5-80% relatief (niet-condenserend)

## 3.7 Emissie en immuniteit

Geluidsemissie

Transportvoorw

Opslagvoorwaa

aarden

rden

< 60 dBA (geluidsdruk), gemeten op een afstand van 1 m van het instrument. Het geluidsniveau kan gedurende korte tijd hoger zijn dan 78 dB terwijl het proces wordt uitgevoerd.

**EMC** De Fluent voldoet aan de eisen voor emissie en immuniteit zoals beschreven in IEC 61326-1 en IEC 61326-2-6. Voor gebruik van de Fluent moet de elektromagnetische omgeving echter worden geëvalueerd. De operator is ervoor verantwoordelijk dat een compatibele elektromagnetische omgeving voor de Fluent kan worden gehandhaafd, zodat de Fluent kan functioneren zoals beoogd. De Fluent is geclassificeerd als APPARATUUR VAN GROEP 1, KLASSE B (CISPR 11).

bij 30 °C (86 °F) of lager



Deze apparatuur is ontworpen voor gebruik in een STANDAARD ELEKTROMAGNETISCHE OMGEVING (IEC 61326-1) en een PROFESSIONELE GEZONDHEIDSZORGINSTELLING (IEC 61326-2-6). De werking is waarschijnlijk onjuist als deze wordt gebruikt in een INDUSTRIËLE ELEKTROMAGNETISCHE OMGEVING (IEC 61326-1) en een THUISZORGOMGEVING (IEC 61326-2-6). Indien het vermoeden bestaat dat de werking wordt beïnvloed door elektromagnetische interferentie, kan de juiste werking worden hersteld door de afstand tussen de apparatuur en de bron van de interferentie te vergroten.

Gebruik de Fluent niet in de nabijheid van bronnen met sterke elektromagnetische straling (bijv. niet-afgeschermde, conventionele RF-bronnen), omdat deze de juiste werking kunnen verstoren.

- **FCC 15** Deze apparatuur is getest en in overeenstemming bevonden met de grenzen voor een digitaal apparaat van klasse B, overeenkomstig deel 15 van de FCCregelgeving. Deze limieten zijn opgesteld om een redelijke bescherming te bieden tegen schadelijke interferentie in een huishoudelijke installatie. Deze apparatuur genereert, gebruikt en kan radiofrequentie-energie uitstralen en kan, indien niet gemonteerd en gebruikt in overeenstemming met de instructies, schadelijke interferentie veroorzaken voor de radiocommunicatie. Er is echter geen garantie dat er geen interferentie zal optreden in een bepaalde installatie. Als deze apparatuur schadelijke interferentie veroorzaakt aan radio- of televisieontvangst, wat kan worden vastgesteld door de apparatuur uit en weer in te schakelen, wordt de gebruiker aangeraden de interferentie te verhelpen door een of meer van de volgende maatregelen te treffen:
  - Draai of verplaats de ontvangstantenne.
  - Vergroot de afstand tussen de apparatuur en de ontvanger.
  - Sluit de apparatuur aan op een stopcontact van een ander circuit dan waarop de ontvanger is aangesloten.
  - Raadpleeg de dealer of een ervaren radio-/tv-technicus voor hulp.



# 4 Beschrijving van de functie

Dit hoofdstuk behandelt het basisprincipe van de Fluent, toont hoe deze is gestructureerd en bevat een functiebeschrijving van de onderdelen.

# 4.1 Overzicht

De Fluent wordt gebruikt voor pipetteertaken met robotarmen. De robotarmen kunnen opzuigen uit en afgeven in verschillende reservoirs, bijvoorbeeld buizen of multititerplaten.

De Fluent is beschikbaar in drie verschillende afmetingen:

- Fluent 480
- Fluent 780
- Fluent 1080



Afb. 17: Instrumentoverzicht (instrument kan van afbeelding afwijken)

A Dek

B Robotarmen

C Opties en apparaten



Een aardbevingsbescherming voor gebieden waar gevaar voor aardbevingen bestaat.

Voor meer informatie, zie paragraaf "Klantsupport" [> 203].

## 4.2 Dek

SegmentenHet Fluent-dek dat de monsterzone van het instrument is, bestaat uit segmenten.<br/>Deksegmenten zijn uitwisselbare dekcomponenten die verschillende afmetingen<br/>en eigenschappen kunnen hebben. VOORZICHTIG! Gebruik het systeem niet<br/>zonder deksegmenten.

# RasternummerDe segmentbreedte wordt uitgedrukt in rasternummers. Een raster is 25 mm breed<br/>en komt overeen met de afstand tussen de positioneerpennen van een segment.



Rasternummers worden ook gebruikt voor het aangeven van segmenten of runners op het dek.

#### 4.2.1 Dragers



Afb. 18: Fluent-dek

А

Runner B Segment

Dragers zijn dekcomponenten voor het vasthouden van laboratoriummateriaal of verbruiksmiddelen op het dek.

Runners zijn dragers die op en van de rastersegmenten schuiven en meestal monsterbuizen of reagensbakken vasthouden.

Segmenten zijn statische elementen die op het dek zijn vergrendeld. Sommige segmenten hebben nesten (nestsegmenten) die laboratoriummateriaal vasthouden, bijv. multititerplaten, deep well-platen of verbruiksmiddelen zoals DiTiboxen. Sommige segmenten hebben rasterpennen (rastersegmenten) voor het inen uitladen van runners.



#### 4.2.2 Dektrays



#### Afb. 19: Dektray

Dektrays, die zich onder de dynamische deksegmenten bevinden, vangen gemorste vloeistof op in het handmatige deklaadbereik. Het systeem moet worden gebruikt met zo veel mogelijk onder het dek gemonteerde dektrays, om alle gemorste vloeistoffen op te kunnen vangen. **VOORZICHTIG! Bedien het systeem niet zonder dektrays en deksegmenten.** 

Uitsparingen in dektrays voor gereedschappen en instrumenten zijn alleen in de kabinetversie toegestaan.

De verhoogde gebieden in elke dektray zijn ontworpen om uitsparingen mogelijk te maken zonder het opvangvolume te beïnvloeden. Deze uitsparingen zijn geschikt voor een afvalkoker door het dek of andere integraties van apparaten bij gebruik van de op een kast gemonteerde versie. Voor toegangspunten door het dek kunnen trays worden geplaatst op een plank in de kast onder geïntegreerde apparaten.



Afb. 20: Dektrays onder de deksegmenten

Dektrays zijn niet aanwezig waar de RGA toegang tot een apparaat onder het dek vereist. Bij het instrument wordt een set dektrays geleverd. De dektrays kunnen worden gewassen of vervangen indien nodig. Zie paragraaf "Einde van de dag" [> 119].



## 4.2.3 Plaatsing op segment



Afb. 21: Vergrendelpennen en positioneerpennen

A Vergrendelpennen B Positioneerpennen

De Fluent gebruikt pennen om runners, adapters of opties correct op een segment te positioneren. De runners zijn ontworpen om op de pennen te schuiven. Het positioneren kan dan worden gecontroleerd door het lezen van het rasternummer aan de voorzijde van het instrument. De vergrendelpennen houden de runner in hun positie.



#### 4.2.4 Segmentpositie



Afb. 22: Zijposities en rasterposities

Zijposities **B** Rasterpositie

De zijposities (AB, YZ) kunnen worden gebruikt om laboratoriummateriaal door de RGA te plaatsen.



Met de FCA of MCA is pipetteren in zijposities niet mogelijk.

De genummerde rasterposities (1–n) zijn toegankelijk voor pipetteerarmen. Bij configuraties met meerdere armen zijn echter niet alle numerieke rasters voor alle pipetteerarmen toegankelijk. Beperkingen kunnen afhankelijk van de instrumentarmconfiguratie van toepassing zijn.

# 4.3 Robotarmen

А

Fluent kan worden uitgerust met meerdere robotarmen:

- Flexibele kanaalarm (FCA)
- Multikanaalarm (MCA 96 en MCA 384)
- Robotgrijperarm (RGA)

De robotarmen kunnen worden uitgerust met meerdere armaccessoires.



### 4.3.1 Flexibele kanaalarm (FCA)



Afb. 23: Flexibele kanaalarm



Als geen residu-overdracht is toegestaan, wordt het gebruik van wegwerppunten met filters sterk aanbevolen.

De FCA (A) is uitgerust met pipetteerpunten en kan vloeistofverwerking voor tot 8 aparte kanalen besturen.

De FCA die is geconfigureerd met DiTi-adapters heeft een optionele FCA-grijper die bepaalde bewegingen mogelijk maakt—zie "FCA-grijper" [▶ 61].

#### 4.3.1.1 FCA met vloeistofsysteem (Liquid FCA)

De FCA met een vloeistofverplaatsingssysteem is gevuld met systeemvloeistof die wordt verzorgd door spuitpompen. Deze wordt gebruikt voor het pipetteren van vloeistoffen met verschillende volumebereiken afhankelijk van de gebruikte punten en de injectiespuiten. De Liquid FCA kan worden geconfigureerd met verschillende vaste punten, wasbare punten of wegwerppunt-adapters.



Tecan adviseert het gebruik van gede-ioniseerd water als systeemvloeistof.

#### 4.3.1.2 FCA met luchtsysteem (Air FCA)

De FCA met een air-displacementsysteem wordt gebruikt voor pipetteren van vloeistoffen door een zuiger in het pipetteerkanaal te bewegen. De Air FCA wordt geconfigureerd met wegwerppunt-adapters.



#### 4.3.2 Multikanaalarm 384 (MCA 384)



Afb. 24: Multikanaalarm



Als geen residu-overdracht is toegestaan, wordt het gebruik van wegwerppunten met filters sterk aanbevolen.

De MCA (A) is een robotarm met een multikanaal-pipetteerkop. Het opzuigen en afgeven gebeurt bij alle pipetteerkopkanalen tegelijkertijd. De kopadapters van de pipetteerkoppen kunnen worden vervangen. Met de verschillende typen kopadapters zijn verschillende pipetteerformaten mogelijk:

- Compatibel met MCA DiTi-portfolio
- Pipetteren tot 125 µl
- Pipetteerbereik tot 500 µl met optionele/uitgebreide volumeadapter
- Geen vloeistofniveaudetectie
- MCA384 met 384 wegwerpnaalden
- MCA384 met 96 wegwerpnaalden (adapterplaat)
- MCA 384 met 384 vaste, wasbare naalden
- MCA 384 met 96 vaste, wasbare naalden

#### 4.3.3 Multikanaalarm 96 (MCA 96)

De MCA 96 is een robotarm met een multikanaal-pipetteerkop. Het opzuigen en afgeven gebeurt bij alle 96 pipetteerkopkanalen tegelijkertijd. De MCA 96-arm heeft de volgende belangrijke kenmerken:





Afb. 25: MCA 96-kop/grijper

- Compatibel met geleidende FCA-DiTi-portfolio tot en met 1000 μl
- Pipetteerbereik van 1 µl tot 1000 µl
- · Voer een (capacitieve) vloeistofniveaudetectie (cLLD) uit
- Voer een gedeeltelijke DiTi-opname en offset-pipettering uit (bijv. voor verdunning) met behulp van naaldproducten met de Tecan-combitray)
- Optionele grijper voor eenvoudige taken met betrekking tot de verwerking van laboratoriummateriaal

*De MCA 96-kop werkt met geleidende wegwerpnaalden op de MCA 96 voor compatibiliteit met cLLD.* 

De cLLD op de MCA 96 is alleen geverifieerd voor gebruik in bakken.

De eerste naald die een vloeistof aanraakt, activeert het cLLD-signaal. Ongelijke vulhoogten (bijv. in handmatig gevulde wells van een multititerplaat) kunnen leiden tot onbedoelde verwerking en foutieve resultaten.

6

Gebruik alleen cLLD-compatibele dragers en cLLD-compatibel laboratoriummateriaal in combinatie met de MCA 96. Gebruik compatibele trays voor gedeeltelijk oppakken van DiTi's: Tecan SLAS-tray Raadpleeg de referentiehandleiding (Ref. [4]) voor meer informatie.

De optionele, in het veld bij te werken grijper op de MCA 96 is geschikt voor eenvoudige taken met betrekking tot de verwerking van laboratoriummateriaal.



De grijper kan bijvoorbeeld worden gebruikt voor de volgende taken:

- Multititerplaten en DWP's in horizontale positie verplaatsen naar doelposities op de werktafel
- Werken met gestapelde DiTi-trays door lege lagen weg te gooien
- Multititerplaatdeksels plaatsen

De MCA-grijper is gebouwd als een submodule op de MCA-kop, dus het is een compact, kostenefficiënt alternatief voor een speciale robotgrijperarm (RGA). De MCA 96-grijper kan geen complexere taken voor de verwerking van laboratoriummateriaal uitvoeren, zoals:

- · Toegang tot hotels
- · Werken met laboratoriummateriaal in verticale positie
- Taken met rotaties

Voor deze complexe taken moet nog steeds een speciale robotgrijperarm (RGA) aan het systeem worden gekoppeld. De MCA 96-grijper bestaat uit hetzelfde vingerwisselsysteem als de RGA-grijper.

#### 4.3.4 Robotgrijperarm (RGA)



Afb. 26: Robotgrijperarm

De RGA (A) is een robotarm met een grijperkop en grijpervingers. De RGA transporteert multititerplaten en ander laboratoriummateriaal tussen dekposities, randapparaten en de opslag voor laboratoriummateriaal:

Een "Robotic Gripper Arm" (RGA standaard Z) met standaard hoogte heeft toegang tot objecten op het dek of op het lagere dek.

Een "Robotic Gripper Arm" (RGA standaard Z) heeft toegang tot objecten op het dek, op het lagere dek en eronder.

#### 4.3.4.1 Robotgrijperkop

De RGA kan worden uitgerust met twee verschillende, optionele robotgrijperkoppen.

De normale grijperkop biedt de keuze uit grijpervingers die handmatig kunnen worden uitgewisseld.



Het "Finger Exchange System" (FES) biedt automatische vingeruitwisseling met een grote keuze aan grijpervingersets. Vingersets zijn bevestigd op een dokstation, gemonteerd op een standaard nestsegment. Vingersets worden door de robotarm automatisch opgenomen en geplaatst. Vingertype en vingerwissel worden bewaakt. Sommige of alle vingers kunnen worden gebruikt binnen een enkele methode.

#### 4.3.5 Armaccessoires



#### Afb. 27: Armaccessoires

A Vaste punten

**C** Grijpervingers

- B Wegwerppunten
  - Barcodescanner voor de RGA

#### 4.3.5.1 Vaste punten



Als geen residu-overdracht is toegestaan, wordt het gebruik van wegwerppunten met filters sterk aanbevolen.

D

Wasbare, herbruikbare punten voor opzuigen en afgeven zijn beschikbaar voor de FCA en MCA.

#### 4.3.5.2 Wegwerppunten

Punten zijn beschikbaar in trays of boxen (enkel of genest) afhankelijk van het type. Punten worden na opzuiging afgevoerd of opnieuw in rekken aangebracht. Punten worden met het wegwerppunt-uitwerpsysteem afgevoerd in een afvoergoot gemonteerd op een deksegment.

#### 4.3.5.3 Grijpervingers

Er zijn verschillende typen grijpervingers beschikbaar voor de normale grijperkop, RGA, en de MCA 96-grijper, beide voorzien van het vingerwisselsysteem (FES).

**Excentrische** grijpervingers Excentrische grijpervingers transporteren multititerplaatgebaseerde objecten binnen en buiten het pipetteerbereik. Plaatobjecten van opzij pakken. Er zijn twee varianten beschikbaar:

- Vingers van standaardlengte voor het laden van multititerplaten in hotels en apparaten.
- Excentrische, lange vingers voor het laden van diepere apparaten zoals de bewaakte incubator met 4-sleuven en celplaat.



MCA 96 excentrische grijpervingers	De excentrische grijpervingers van de MCA 96 transporteren multititerplaatgebaseerde objecten binnen het pipetteerbereik. Dit vingertype kan plaatobjecten onder de grijperkop van de MCA 96 vastgrijpen. (Dit vingertype is NIET compatibel met de RGA-grijper.)
Centrische grijpervingers	Centrische grijpervingers transporteren multititerplaatgebaseerde objecten binnen en onder het pipetteerbereik. Plaatobjecten van boven pakken. Alleen van toepassing op de normale grijperkop, RGA.
Buisvingers	Buisvingers transporteren buisgebaseerde objecten binnen en onder het pipetteerbereik. Alleen van toepassing op de normale grijperkop, RGA.

#### 4.3.5.4 Barcodescanner

De RGA kan worden uitgerust met een horizontale scanner voor barcodes of multititerplaten en DiTi-boxen.



De veiligheidsinstructies van de laserklasse moeten zorgvuldig worden gelezen en opgevolgd.

Zie ook de handleiding van de fabrikant van de barcodescanner.

# 4.4 Vloeistofsysteem (Liquid FCA)



Afb. 28: Vloeistofsysteem (Liquid FCA)

Het vloeistofsysteem is ontworpen voor het efficiënt wassen, in- en uitwendig, van vaste pipetteerpunten.





# 4.5 Wassysteem (MCA 384)



Afb. 29: Wassysteem (MCA)

Het wasblok (A), gemonteerd op het MCA 384-segment, dient om de naalden van de vastenaaldadapter na elke pipetteercyclus te wassen.

# 4.6 Opties en apparaten



Voorbeeld van een lijst met opties en apparaten van derden voor de integratie op Fluent. Bepaalde opties van Tecan en apparaten van derden die met de Fluent kunnen worden gebruikt zijn uitsluitend voor onderzoekstaken bedoeld ('researchuse-only', RUO).

*In deze paragraaf zijn opties en apparaten voor alleen-onderzoek gemarkeerd met een asterisk (\*).* 

Voor meer informatie, zie paragraaf "Beoogd gebruik" [> 8].

**Passieve opties** 

- Hotel (plaatopslag)
- Kabinet
- Stofafdekking
- FCA-grijper

**Actieve opties** 

- HEPA-kapFluent Stacker
- MIO2
- Te-Shake
- Te-VacS
- Fluent-carrousel
- Was- en navulcentrum (WRC) torens (bijv. MCA-wasstation)
- Resolvex i300
- Doorpriknaalden en buisrotator
- FRIDA Reader

Voor meer informatie, zie paragraaf "Referentiedocumenten" [> 9].

• Waseenheden gebaseerd op HydroControl



- Balansen gebaseerd op de MT-SICS niveau 1 standaard
- SiLA-compatibele apparaten\*
- Agilent-sealer\*
- Inheco ODTC
- Inheco Heating Cooling met MTC/STC-controller
- Cytomat 10\*, 20\*, 200\* en 6000\*

#### Barcodelezers

Fluent ID-buisbarcodescannerBarcodelezers van de Keynence-serie BL-1300

Lezers

- Tecan-lezers bestuurd door Magellan
- Spark en SparkControl Magellan\*
- Ziath 2D flatbed-lezer\*



Zie ook de handleidingen van de fabrikant van de optie, het apparaat of het externe apparaat. De instructies moeten zorgvuldig worden gelezen en opgevolgd.

#### 4.6.1 Fluent ID-buisbarcodescanner



#### Afb. 30: Fluent-ID

- A Scannerbehuizing
- C Laadbereik
- E Reflector

- B Laserbarcodescanner
- D Leds

De Fluent ID is een optionele module die kan worden geïntegreerd om barcodelabels te scannen, als buisrunners op het dek worden geladen. Elke Fluent ID-module bevat zes specifieke rasterposities voor het laden en scannen van de barcodelabels van maximaal zes runners. De reflector wordt gebruikt voor het detecteren van lege buisposities in een runner. Met een grafische interface op het touchscreen kan door de Fluent ID-bediening worden genavigeerd.

De laserstraling van de barcodescanner is een laagvermogen, gecollimeerde bundel in het zichtbare spectrum met de volgende eigenschappen:



- Golflengte: 655 nm
- Impulsduur: 150 µs
- Maximaal vermogen van energie-output: 1,0 mW

#### 4.6.1.1 Met Fluent ID compatibele buisrunners

De Fluent ID-buisrunners zijn steeds voor een bepaald buistype ontworpen:

- Runner met 32 posities voor buizen met 10 mm diameter
- Runner met 32 posities voor buizen met 13 mm diameter
- · Runner met 26 posities voor buizen met 16 mm diameter
- Runner met 32 posities voor 2ml Eppendorf Safe-Lock-buizen



Optionele pluggen kunnen worden gebruikt voor het blokkeren van twee posities van een 26-positierunner om deze te gebruiken als een 24-positierunner waardoor parallel pipetteren uit buizen in veelvouden van acht mogelijk is.

#### 4.6.2 FCA-grijper

#### **Overzicht**



Afb. 31: FCA-grijper

De FCA-grijper is een optie voor de FCA geconfigureerd met DiTi-adapters waarmee — naast pipetteren — de FCA nog andere bewegingen van laboratoriummateriaal kan maken. Tijdens de cyclus kan de FCA de FCAgrijpervingers automatisch oppakken en neerlaten. **4 - Beschrijving van de functie** Opties en apparaten



FCAgrijpervingers



Afb. 32: FCA-grijpervingers

De FCA-grijpervingers moeten na 2 jaar of 20000 gebruikscycli worden vervangen (een cyclus gedefinieerd als oppakken, gebruiken en parkeren). De cycli worden bewaakt met een teller zoals gedefinieerd in de Fluent Control-software.



Afb. 33: FCA-grijper dokstation-nest

Het FCA-grijper dokstation-nest wordt gebruikt voor het opbergen van FCAgrijpervingers. Het kan als een standaard multititerplaatnest op een deksegment worden bevestigd.

FCA-grijper dokstation-nest



#### 4.6.3 Mix & Pierce

Het Mix & Pierce-werkstation van de Fluent is ontworpen voor toepassingen die middels doorprikken vloeistof van en naar met rubberdop afgesloten monsterbuizen overdragen zonder dat de rubberdoppen hoeven te worden verwijderd.



De septa van de buis kan slechts eenmaal worden doorgeprikt. Herhaaldelijk doorprikken van dezelfde buis wordt niet ondersteund.

6



BD Vacutainer®-buisjes met Hemogard-sluiting en Greiner Vacuette®-buisjes met niet-geribbelde trekdop en veiligheidsdraaidop zijn getest op doorprikken. Doorprikken van buizen met rubberen stoppen/pluggen wordt niet ondersteund.

Doorpriknaalden moeten regelmatig worden vervangen. Het gebruik is geverifieerd voor 20.000 doorprikkingen/naald.

Het Mix & Pierce-werkstation van de Fluent is geconfigureerd met tot 2 vloeistof-FCA's, een diep wasstation en tot 4 buisrotators op de Fluent-basiseenheid. Buisrotators kunnen op elke Fluent-basiseenheid worden geïntegreerd en ondersteunen het lezen van barcodes, het mengen, doorprikken en aliquoteren van monsters. Voor meer informatie over ondersteunde buistypen, zie paragraaf "Buisrotator-runners" [> 64].

De verwerking kan in de volgende stappen worden verdeeld:

- 1. Barcode scannen tijdens laden van buizen
- 2. Mengen van buisinhoud
- 3. Doorprikken en vloeistof verwerken met de FCA in de buisrotator met doorpriknaalden
- 4. Wassen en ontsmetten van doorpriknaalden in het diepe wasstation en ontsmettingsbakken
- 5. Herhaling van stap 2 en voortzetting
- 6. Aan het einde van het script wordt een uitgebreide spoeling van het wasstation aanbevolen, met inbegrip van de centrale afvalreinigers voor en achter.

#### 4.6.3.1 Buisrotator

Het hoofddoel van de buisrotator is om de vloeistofinhoud van de buizen te mengen en te dienen als drager voor doorprik- en pipetteeracties. Een enkele buisrotator heeft een capaciteit van 5 buisrotator-runners met elk 24 buizen (bijv. 120 buizen totale capaciteit).

Het apparaat bevat de volgende subcomponenten:

- Een geïntegreerde buisbarcodescanner voor het scannen van de monsterbarcodes tijdens laden
- Een optioneel diep wasstation met diepe bakken voor het ontsmetten van doorprikpunten en een foutbuisdrager. De houder voor foute buizen kan worden gebruikt voor het bewaren van monsters bij doorprikfouten. Het wasstation wordt naast de rotatietrommel geplaatst.



- Een roterende trommel met buisneerhouder die plaats biedt aan maximaal vijf buisrotator-runners. De trommel voert het mengen van monsters middels 360°rotatie of oscillatie onder verschillende hoeken en snelheden uit. De neerhouder (afdekking) ondersteunt het doorprikproces.
- De buisrotator met buisrotator-runners ondersteunt capacitieve vloeistofniveaudetectie voor en na de opzuiging en na vloeistofafgifte (vloeistofaankomstcontrole) door gesloten buizen (optionele instelling).
- De buisrotator is geïnstalleerd door de technicus en mag niet door de keyoperator of gebruiker worden verplaatst.

#### 4.6.3.2 Buisrotator-runners

Buisrotator-runners zijn ontworpen voor het gebruik op de buisrotator-runner en ter ondersteuning van het doorprikken. Er bestaan verschillende buisrunners om het doorprikken voor verschillende buistypen mogelijk te maken:

- 13 x 75 mm BD buisrotator, 24 buisposities
- 13 x 100 mm BD buisrotator, 24 buisposities
- 13 x 75 mm Greiner buisrotator-runner, 24 buisposities
- 13 x 100 mm Greiner buisrotator-runner, 24 buisposities
- 16 x 100 mm buisrotator-runner, 24 buisposities



*Meervoudige afgifte wordt alleen ondersteund voor Greiner Vacuette*®*-buisjes met een niet-geribbelde trekdop en veiligheidsdraaidop.* 

Productlijn	Buis			Runner	
	Diame- ter [mm]	Lengte [mm]	Definitie laboratorium- materiaal	Compatibele runner	Brug- kleur
Vacuette van Greiner	13	100	13 x 100 mm Greiner Vacuette met septum	1x24 13 x 100 mm Greiner buisrotator-runner	grijs
	13	75	13 x 75 mm Greiner Vacuette met septum	1x24 13 x 75 mm Greiner buisrotator-runner	
	16	100	16 x 100 mm Greiner Vacuette met septum	1x24 16 x 100 mm buisro- tator-runner	zwart
BD Vacutainer	13	100	13 x 100 mm BD Vacu- tainer met septum	1x24 13 x 100 mm buisro- tator-runner	wit
	13	75	13 x 75 mm BD Vacu- tainer met septum	1x24 13 x 75 mm buisrota- tor-runner	
	16	100	16 x 100 mm BD Vacu- tainer met septum	1x24 16 x 100 mm buisro- tator-runner	zwart

#### Tab. 5: Buis- en runnercompatibiliteit





#### 4.6.3.3 Doorprikpuntbescherming



Afb. 34: Doorprikpuntbescherming

De doorprikpuntbescherming is een dop die wordt gebruikt voor het afdekken van de scherpe punt van de piercingpunten tijdens het vervangen van de punt en het verhelpen van storingen. Hij beschermt de gebruiker tegen letsel en de punten tegen beschadiging.



De doorprikpuntbescherming is bedoeld voor eenmalig gebruik. Na gebruik moeten alle doorprikpuntbeschermingen worden afgevoerd in het biologische afvoerreservoir.



#### 4.6.3.4 Doorprikpunt-verwijderingsgereedschap

Afb. 35: Doorprikpunt-verwijderingsgereedschap

Het doorprikpunt-verwijderingsgereedschap wordt gebruikt voor het terugtrekken van een doorprikpunt die vast zit in een buis en niet meer via de software kan worden teruggetrokken.

#### 4.6.4 Frida Reader



Afb. 36: Frida Reader



De Frida Reader is bedoeld voor de geautomatiseerde kwanticering en normalisering van nucleïnezuren. Monsters die met de Frida Reader worden gemeten, moeten bij 4°C worden gekoeld, anders kan verdamping van het monster de meetresultaten beïnvloeden.

# **A VOORZICHTIG**

#### Trillingen kunnen leiden tot onjuiste resultaten!

Trillingen van de monsterdruppel kunnen leiden tot onjuiste meetresultaten en afbreuk doen aan de veiligheid of de klinische toestand van het patiëntenmonster.

- Een stabiele ondergrond is een eerste vereiste voor een geschikte opstelplaats.
- Tijdens metingen van de Frida Reader zijn geen interne of externe trillingsbronnen in de buurt toegestaan.
- Vermijd bronnen met resonantiefrequentie. Met name trillingen rond 36 Hz (2160 omw/min) en rond 42 Hz (2520 omw/min) moeten worden vermeden, omdat dit resonantiefrequenties zijn van een hangende druppel.

# 

#### Ruimteverlichtingen kan leiden tot onjuiste meetresultaten!

Ruimteverlichting boven de module kan de meting storen kan leiden tot onjuiste meetresultaten en afbreuk doen aan de veiligheid of de klinische toestand van het patiëntenmonster.

 Het robotsysteem moet zijn voorzien van een niet-transparante afdekking, een voor- en achterpaneel, om omgevingslicht op de meetpositie van de Frida Reader te voorkomen.

#### 4.6.5 Fasescheider

De fasescheider is ontworpen om scheidingsfasen te detecteren tussen vloeistoffen met verschillende viscositeit. Als zodanig is deze onafhankelijk van de scheidingsfase die zichtbaar is van buiten het laboratoriummateriaal. De fasescheider kan worden gebruikt in toepassingen die de probleemloze overdracht van een vloeibare fase van een bron naar doellaboratoriummateriaal vereisen.

Hieronder wordt een standaardwerkwijze weergegeven:

- ✓ Centrifugeren van het bronlaboratoriummateriaal met een vloeibaar mengsel om een afzonderlijke fase tussen vloeistoffen te genereren.
- De vloeistoffen moeten verschillen in viscositeit om fasevorming tijdens het centrifugeren mogelijk te maken.
- Plaats de buisjes zonder doppen op het Fluent-platform en gebruik een barcodescanner (bijv. Fluent ID) voor volledige traceerbaarheid. Zorg ervoor dat de laag tussen de fasen/vloeistoffracties tijdens het laden niet wordt verstoord.
- Start het protocol dat is gedefinieerd voor scheiding van de fracties. De fasescheidingsfunctie van de Air FCA detecteert de fase tussen de vloeistoffen en start de overdracht van de gewenste fractie naar het doellaboratoriummateriaal.



Er kan meer dan één fase uit het bronlaboratoriummateriaal worden geëxtraheerd. Het verwijderen van ten minste een deel van de bovenste fase is vereist om te voorkomen dat er vloeistof op het platform wordt gemorst (overstromen van laboratoriummateriaal) en om mogelijke verontreiniging van het pipetteerkanaal boven de wegwerpnaald tijdens fasedetectie te voorkomen. Baadpleeg de referentiebandleiding voor technische specificaties. Baadpleeg de

Raadpleeg de referentiehandleiding voor technische specificaties. Raadpleeg de softwarehandleiding van de toepassing voor meer informatie over . Zie "Referentiedocumenten" [> 9].

De vloeibare fasen moeten duidelijk worden gescheiden. Voor de scheiding van volbloed in plasma en bloedcellen is de kwaliteit en voorbehandeling van de monsters van essentieel belang. Parameters die van invloed kunnen zijn op de fasedetectie in bloedmonsters zijn de kwaliteit van het monster (lipemie, hemolyse), de opslagduur, de opslagtemperatuur, de transportomstandigheden, centrifugeomstandigheden (tijd, rcf, temperatuur, helling, rotortype), vervorming van de fase na centrifugatie enz.

Voor optimale resultaten van fasescheiding moeten bloedmonsters zo snel mogelijk na afname worden verwerkt. De monsterbehandeling en opslagomstandigheden moeten voldoen aan de aanbevelingen van de fabrikant van de desbetreffende buisjes.

Centrifugeren bij 2500 rcf gedurende 10 minuten bij kamertemperatuur met een vertraagde helling leidt tot een probleemloze fasescheiding voor plasmamonsters (interne testomstandigheden van Tecan).

Om te voorkomen dat er tijdens het opzuigen en de fasedetectie vloeistof uit de buisjes loopt, mogen de buisjes niet tot de rand worden gevuld. Voor fasedetectie is gewoonlijk een snelle neerwaartse beweging in het buisje vereist in combinatie met een langzame opzuigsnelheid, wat leidt tot een stijging van het vloeistofniveau tijdens detectie.

Voor buisjes met een hoog vulvolume wordt aanbevolen om vloeistof van de bovenkant te verwijderen voordat de fasedetectie wordt gestart.



#### 4.6.6 Resolvex i300



Afb. 37: Resolvex i300-werktafelmodule

De Resolvex i300 is een module die in het -platform kan worden geïntegreerd. Het -platform kan vervolgens vloeistoffen verwerken en laboratoriummateriaal rechtstreeks vanaf een werktafel overbrengen naar de Resolvex i300-module binnen hetzelfde softwareframework.

De Resolvex i300 is een laboratoriuminstrument waarmee een reeks laboratoriumprocessen kan worden geautomatiseerd met behulp van de volgende hoofdfuncties:

- Filterkolommen (volledig of gedeeltelijk geladen) onder druk zetten met lucht of stikstof
- Stapelen en ontstapelen van filterplaten op verzamelplaten met een geïntegreerde lift voor laboratoriummateriaal
- Bescherming van monsters tegen kruisbesmetting tijdens het stapelen en ontstapelen van filter- en verzamellaboratoriummateriaal met behulp van een geïntegreerde druppelbescherming
- Aftappen van vloeistoffen in het filter of verzamelen van laboratoriummateriaal (optioneel)
- Verdamping van vloeistoffen uit het laboratoriummateriaal met verwarmd gas, zoals lucht of stikstof (optioneel)



Gebruikelijke toepassingsworkflows van de Resolvex i300 omvatten de volgende laboratoriumprocessen:

- SPE-processen (Solid phase extraction, vaste-fasextractie) voor workflows voor monstervoorbereiding voor massaspectrometrie
- Ondersteunde vloeistofextractie (Supported liquid extraction, SLE)
- Andere filtratie-, zuiverings- of concentratieworkflows met positieve druk

Raadpleeg voor meer informatie en voordat u werkzaamheden uitvoert met de Resolvex i300 de bedieningshandleiding voor de Resolvex i300. Zie "Referentiedocumenten" [> 9].



5 - Besturingselementen Bedieningselementen

# **5** Besturingselementen

# 5.1 Bedieningselementen



Het touchscreen toont methoden en beschrijvingen waarmee de operator het instrument kan bedienen.



# 5.2 Gebruikersinterface



C Display/optie/actietoetsen

Door de gebruikersinterface van FluentControl heeft de operator toegang tot de methodecycli voor bediening en systeemverzorging.

#### 5.2.1 Navigatiepad

Gebruik het navigatiepad om de hiërarchie van FluentControl te begrijpen en erin te navigeren.

Tab. 6: Navigatiepadtoetsen

Toets	Naam	Functie
۵	Home	Indrukken om terug te keren naar de startpagi- na.
System Care 🛞	Navigatiepaneel	Voor het weergeven van actuele en eerdere se- lecties.
<	Menu-uitbreiding	Druk op de menu-uit- breidingstoets voor toe- gang tot opties zoals lichtbediening en om de operators te wisselen.


### 5.2.2 Werkbereik

Toegangsmethoden en -beschrijvingen via het werkbereik van de gebruikersinterface. Details over de status van de methodecyclus worden ook hier weergegeven.

Tab. 7: Werkbereik-toetsen

Toets	Naam	Functie
	Cyclus	Indrukken om de gese- lecteerde methode te starten.
+	Toevoegen	Indrukken om meer methoden aan uw snel- startlijst toe te voegen.
Assay 1 unknown	Geselecteerde methode	De actueel geselecteer- de methode die wordt uitgevoerd, als Volgen- de wordt ingedrukt.
Assay 3 unknown	Beschikbare methode	Een methode die kan worden geselecteerd door erop te klikken.
Start of Day Due: is not executed	Snelstart-toets	Indrukken om de gese- lecteerde methode on- middellijk te starten.

Tab. 8: Werkbereikdisplay

Display	Displayfunctie
00:16:52 10:09 10:43 Tot Time	Toont de status en de resterende tijd voor de methodecyclus.
Assay 1 is ready to be started.	Beschrijving van de actueel geselec- teerde methode of aanvullende infor- matie over de actuele actie.

### 5.2.3 Display, optie en actietoetsen

Tab. 9: Display, optie en actietoetsen

Toets	Naam	Functie
✓ Ok	Ok	Indrukken om te bevestigen.



Toets	Naam	Functie
<b>X</b> Cancel	Annuleren	Indrukken om te annuleren.
Continue	Verder	Indrukken om verder te gaan.
<b>II</b> Pause	Pauze	Indrukken om een pauze aan te vra- gen aan het einde van de actuele actie.
U Stop	Stop	Indrukken om een cyclus onmiddellijk te stopen, zelfs midden in een actie- ve actie. Indien mogelijk, biedt het systeem de mogelijkheid om de cyclus te her- stellen of voort te zetten.
ffffffffffffffffffffffffffffffffffffff	Verwijderen	Indrukken om de methode uit het snelstartaanzicht te verwijderen.
View Mode	Aanzichtmodus	Indrukken om te wisselen tussen lijstaanzicht en snelstartaanzichten.
Sort by	Sorteren naar	Indrukken om in het methodecyclus- display te wisselen tussen alfabeti- sche sortering en sortering naar ac- tualiteit.

### 5.2.4 Methodeherstel-toetsen

Tab. 10: Display, optie en actietoetsen

Toets	Naam	Functie
Discard	Afvoeren	Indrukken om een herstelde metho- destatus af te voeren.
$\rightarrow$	_	Indrukken om naar het volgende scherm te wisselen.



Toets	Naam	Functie
<b>C</b> Recovery Point	Herstelpunt	Indrukken om terug te keren naar het vorige scherm ("Herstelpunt").
Run Recovery	Cyclusherstel	Indrukken om de cyclus voort te zet- ten.

### 5.2.5 DeckCheck-toetsen

Tab. 11: DeckCheck-toetsen

Toets	Naam	Functie
Left	Linker camera	Toont de camera-afbeelding van de linker camera (alleen Fluent 780/1080). Een uitroeptreken ver- schijnt op het pictogram als een af- wijkende dekindeling met deze ca- mera werd geregistreerd.
Center	Middencamera	Toont de camera-afbeelding van de overzichtscamera in het midden. Een uitroeptreken verschijnt op het picto- gram als een afwijkende dekindeling met deze camera werd geregistreerd.
Right	Rechter camera	Toont de camera-afbeelding van de rechter camera (alleen Fluent 780/1080). Een uitroeptreken ver- schijnt op het pictogram als een af- wijkende dekindeling met deze ca- mera werd geregistreerd.
Pause Alternate	Wisselen stoppen	Scherm wisselt tussen referentie- en live-afbeeldingen: Druk op deze toets, terwijl een refe- rentie- of live-afbeelding wordt weergegeven om deze afbeelding statisch te houden.
Resume Alternate	Wisselen voortzetten	Afbeelding is statisch: Druk op deze toets om wisselen tus- sen referentie- en live-afbeeldingen voort te zetten.



Toets	Naam	Functie
Check	Check	Activeert een hercontrole van het systeem — bijvoorbeeld nadat cor- recties zijn uitgevoerd. Er wordt ver- zocht de deur te sluiten. Bij een 3-armig systeem moet de middenarm bewegen: Als de deur niet gesloten is, wordt de controle uitgevoerd, maar blokkeert de middelste arm één camera.
Ignore & Continue	Ignore & Continue	Verschijnt alleen, indien geconfigu- reerd voor die opdracht in de metho- de. Hiermee kunnen gemarkeerde afwijkingen worden genegeerd, terwijl het script verder wordt uitgevoerd.
Continue	Continue	Verschijnt, als alle afwijkingen zijn opgelost of als het systeem geen af- wijkingen heeft gevonden en de optie <b>altijd weergeven</b> werd geselecteerd voor de opdracht. Hierdoor zijn sub- tiele kleurveranderingen mogelijk die het menselijk oog wel ziet, maar het systeem niet herkent.

### Tab. 12: Displays

Display	Beschrijving	Functie
	Referentie-afbeelding	De referentie-afbeelding wordt op- geslagen in de scriptopdracht die de gewenste dekindeling weergeeft.
LIVE	Live-afbeelding	De live-afbeelding van de camera's opgenomen tijdens het uitvoeren van het script.
	Afwijking (verschil met referentie-af- beelding)	Rode vierkanten markeren bereiken op live-afbeeldingen waar afwijkin- gen van de referentie-afbeeldingen werden vastgesteld. Het gemarkeer- de bereik kan meer dan één fout be- vatten.





## 5.3 Foutsignalen en instrumentstatus

Afb. 40: Statuslampen

A Voeding-statuslamp B



De statuslamp geeft de instrumentstatus aan middels gekleurde, continu brandende of knipperende lampen. De bovenste statuslamp is alleen actief, als software actief is.

Bovenste statuslamp

Tab. 13: Lichtsignalen van statuslampen

Sig- naal	Kleur	Modus	Instrumentstatus
	_	uit	Het instrument is uitgeschakeld (los- gekoppeld van de voeding).
	wit	"hartslag"	Het instrument is ingeschakeld (be- sturingssoftware verbonden, modules nog niet geïnitialiseerd).
	wit (alleen voedings- lamp)	continu	Instrumentstatus "voeding aan" (be- sturingssoftware is niet verbonden).



Sig- naal	Kleur	Modus	Instrumentstatus
	kleurschema	"hartslag"	Stationaire modus
	van Fluent- Control ge- bruikersinter-		Alle modules worden geïnitialiseerd; het instrument is klaar voor het uit- voeren van een methode.
	lace		Na ongeveer een uur in stationaire modus, schakelt het instrument naar stand-bymodus.
			Stand-bymodus
			Alle assen worden geremd. De ar- men zijn niet in ZeroG en kunnen niet handmatig worden bewogen. Om het instrument te activeren, een methode uitvoeren of de key-operator vragen om de Move Tool voor ZeroG modus te selecteren.
	geel	continu	Teach-modus
			Het instrument "leert" posities.
			In deze modus kan de gebruiker de robotarmen handmatig bewegen.
	groen	continu	Een methode (script of proces) is actief.
			Dit is de reguliere <b>"productie -mo-</b> dus.
	rood	knipperend	Fouttoestand
			Het scherm van de besturingscom- puter of het touchscreen toont een foutmelding.
	kleur kan	knipperend	Gebruikersprompt
	door de ge- bruiker wor- den geconfi- gureerd		Systeem wacht op een gebruikers- handeling.
	groen	knipperend	Actieve stop
			Dit is een bedoelde pauze die wordt geactiveerd door de runtime-control- ler of door het openen van een vei- ligheidspaneel.
			Het instrument pauzeert zodat de gebruiker handelingen aan het dek kan uitvoeren. De operator kan de methode voortzetten.





### 5.4 Fluent ID-statusleds



Afb. 41: Fluent ID-leds

De Fluent ID-leds geven de volgende statussen weer:

Tab. 14: Fluent ID-leds

Sig- naal	Kleur	Modus	Instrumentstatus
	_	uit	Fluent ID is stationair.
	wit	continu	Fluent ID ingeschakeld (maar nog niet geïnitialiseerd).
••••	blauw of aangepaste kleur	knipperend	Klaar voor het in- en uitladen van runners.
	groen	continu	Barcodes succesvol gescand. Runner onder toezicht. Niet uitladen omdat dit de cyclus onderbreekt.
	rood	knipperend	Fouttoestand Foutmelding en vereiste actie worden op het touchscreen weergegeven.



# 6 Bediening

### 6.1 Veiligheidsinstructies voor dit hoofdstuk

### **⚠ VOORZICHTIG**

### Onjuiste resultaten of verontreiniging van instrument!

Onjuiste resultaten of verontreiniging van het instrument kunnen optreden, als de installatiekwalificatie en bedieningskwalificatie niet werden nagekomen of als de gebruiksprocedures in deze bedieningshandleiding niet worden opgevolgd.

- Protocollen met betrekking tot installatiekwalificatie en gebruikskwalificatie zijn beschikbaar en bekend.
- Methoden en processen, inclusief pipettereerparameters moeten door de keyoperator worden gevalideerd.
- Vloeistofniveaudetectie in combinatie met doorpriktoepassingen voor FCA en Air FCA moeten door de key-operator worden gevalideerd.
- Voor volbloedtoepassingen met doorpriknaalden en buisrotator en geconcentreerde corrosieve ontsmettingsoplossingen wordt aanbevolen beide connectoren van het wasstation (bij het wasstation en bij de vloeistofcontainer van het systeem) elke drie maanden te vervangen om verstopping, slijtage en scheuren van de connectors te voorkomen. Potentieel corrosieve vloeistoffen moeten vóór de drainage worden geneutraliseerd of het wasstation moet worden gespoeld met neutrale vloeistoffen, zoals water.
- De operator moet worden geschoold over de bedieningsprocedures, methoden en processen.



### **▲ VOORZICHTIG**

### Biologische en chemische verontreiniging van de gebruiker!

Beschadigde FCA-grijpervingers kunnen platen verliezen. Gevallen platen kunnen door gevaarlijke stoffen verontreiniging veroorzaken.

• Controleer de FCA-grijpervingers na een botsing.

### **⚠ VOORZICHTIG**

#### Scherpe randen en punten!

De doorpriknaalden van het Fluent Mix & Pierce-werkstation hebben scherpe punten en randen die letsel kunnen veroorzaken.

- Beweeg de FCA tijdens het laden van het instrument middels een softwareopdracht naar een veilige positie.
- Dek de doorpriknaalden na een fout af met doorprikpuntbeschermingen, en beweeg de FCA handmatig naar een veilige positie. Zie paragraaf "Doorprikpuntbescherming" [> 65].

### **▲ VOORZICHTIG**

#### Biologische besmetting van het systeem!

In het Fluent Mix & Pierce-werkstation kan bloed de doppen van de buizen besmetten.

- Behandel de buizen voorzichtig.
- Draag beschermingsmiddelen.

### LET OP

#### Storing door corrosieve vloeistoffen!

Drainage van corrosieve geconcentreerde vloeistoffen, zoals 2% bleekmiddel, via het wasstation en de slangen kan storingen in de connectors van het wasstation veroorzaken.

 Potentieel corrosieve vloeistoffen moeten vóór de drainage worden geneutraliseerd of het wasstation moet worden gespoeld met neutrale vloeistoffen, zoals water.

### 6.2 Bedrijfsmodi

De Fluent kan in drie verschillende bedrijfsmodi werken:

Operator

#### Routinemodus

- Normale bedrijfsmodus waarin de toepassing wordt of de routinesysteemverzorgingstaken worden uitgevoerd.
- Fluent wordt bewaakt door de FluentControl-softwareruntimecontroller.



Key-operator	Methode definitiemodus
	• Deze modus wordt gebruikt voor het uitvoeren van speciale taken zoals het aanpassen of instellen van de methode.

#### Technicus Servicemodus

- Deze modus wordt gebruikt voor het uitvoeren van speciale taken zoals controles om de bedrijfsgereedheid te garanderen.
- Een Service-certificaat is vereist om deze modus uit te voeren.

### 

#### Arm botst tegen werkbare objecten aan

Vloeiende armen kunnen handmatig worden bewogen. Zorg ervoor dat de handmatige bewegingen van de armen soepel worden uitgevoerd terwijl de arm wordt vastgehouden. Stoot de armen niet tegen vaste voorwerpen, waaronder de mechanische stop voor het bereik van de armen

### 6.3 Inbedrijfstelling

#### 6.3.1 Het instrument inschakelen

Schakel het instrument als volgt in:

1. Schakel de voeding in met voedingsschakelaar (A) aan de achterzijde van de externe stroomvoorziening.





Als het instrument van stroom wordt voorzien, brandt de bedrijfslamp blauw. Zie paragraaf .

Als de statuslamp niet brandt, de pc starten of de key-operator raadplegen.

2. Start de FluentControl-software. Zie paragraaf "FluentControl starten" [> 83].

### 6.3.2 FluentControl starten

- ✓ De bedrijfsprocedures moeten beschikbaar en bekend zijn.
- ✓ Protocollen met betrekking tot installatiekwalificatie en gebruikskwalificatie zijn beschikbaar en bekend.
- ✓ Systeemverzorging is uitgevoerd.
- ✓ Instrument is ingeschakeld.
- Start de software met Start > Alle programma's > Tecan > FluentControl. Na enkele seconden verschijnt het Start scherm.

### 6.3.3 Gebruikersaanmelding



Gebruikersbeheer is alleen beschikbaar met de optie Fluent Gx Assurancesoftware.

Meld u als volg bij FluentControl aan:

- ✓ Fluent Gx Assurance-software is geïnstalleerd.
- ✓ FluentControl is gestart.
- ✓ Gebruikersbeheer is geactiveerd in FluentControl en het proces is voltooid.
- 1. Selecteer het toegewezen gebruikersprofiel (A).



2. Voer het wachtwoord in op het toetsenbord (B).



3. Druk op **OK** (C).



Bij de volgende aanmelding wordt het instrument automatisch geïnitialiseerd.

### 6.3.4 Segmenten plaatsen

Plaats alle segmenten als volgt:

- ✓ Alle segmenten, dragers, opties en apparaten moeten worden geplaatst overeenkomstig de geselecteerde methode.
- ✓ Segmenten zijn gereinigd en in perfecte toestand.
- ✓ Segmenten worden in de desbetreffende rasterpositie geplaatst.
- 1. Sluit bij het aansluiten van een MCA 384 Active Carrier de kabel aan.



2. Laat het segment neer op de achterzijde van het dek.



- 3. Lijn de achterrand uit ten opzichte van de achterkanaalafdekking of de instrumentuitbreiding.
- 4. Laat het voorste deel van het segment (A) neer.



5. Draai de klemhendel van links naar rechts, naar de gesloten positie. Zie paragraaf "Segment controleren" [▶ 196].

### 6.3.5 Segmenten verwijderen

6

Fluent ID-segmenten zijn niet ontworpen om te worden verwijderd! Ze zijn direct met de instrumentelektronica verbonden. De aansluiting op een elektronisch systeem mag alleen worden uitgevoerd door een gekwalificeerde technicus. Raadpleeg de paragraaf Klantsupport.

Verwijder segmenten als volgt:

- ✓ Alle reagentia, monsters, rekken, runners en platen werden van het segment verwijderd.
- ✓ Er is niets op het segment geplaatst.
- 1. Draai de klemhendel (A) van rechts naar links in de geopende positie.





Het segment is ontgrendeld en de gele markering op de klemhendel is zichtbaar.

2. Duw het segment ongeveer 4 mm naar voren.





3. Til het segment (B) aan de voorzijde op.

Ontkoppel de kabel, alvorens een MCA Active Carrier te verwijderen.



4. Bewaar het segment op een schone en droge plek om schade te voorkomen.

### 6.3.6 Standaard runners inladen

### LET OP

#### Beschadiging door verkeerd in- en uitladen

Schade aan runners en pennen.

- Lijn de runner horizontaal uit ten opzichte van het dek.
- Ondersteun het voorste einde van de runner met een hand.
- Zorg er tijdens uitladen voor dat de runner alle pennen vrijmaakt, alvorens de runner op te tillen.



Laad runners als volgt in:

1. Lijn de runner uit ten opzichte van de desbetreffende rasterpositie (A).



- 2. Duw de runner tegen de stoppositie.
- 3. Zorg ervoor dat de runner het segment veilig vergrendelt.

Dit kan worden gevoeld op de laatste paar millimeter voordat de runner de stoppositie raakt.



Laad runners als volgt uit:

- 1. Trek de runner horizontaal ter hoogte van het dek tot deze volledig uit het laadbereik is verwijderd.
- 2. Ondersteun het voorste einde van de runner met een hand.
- 3. Zorg er voor dat de runner alle pennen vrijmaakt, alvorens de runner op te tillen.



### 6.3.7 Dekindeling controleren

Zorg ervoor dat op het dek geïnstalleerde dragers, laboratoriummateriaal en apparaten overeenkomen met de dekindeling die voor deze methode is gedefinieerd.

### LET OP

### Materiële schade!

Onjuist positioneren van segmenten en laboratoriummateriaal op de werktafel kan leiden tot verbrijzelen van de armen.

- Zorg er altijd voor dat de dekconfiguratie en het geladen laboratoriummateriaal overeenkomen met de FluentControl-werktafelconfiguratie.
- Zorg er altijd voor dat het laboratoriummateriaal correct in de nesten is aangebracht. Zie "Laboratoriummateriaal positioneren" [▶ 171].

### LET OP

### Magnetisch veld veroorzaakt storing!

Een sterk magnetisch veld (noordpool omhoog) bij de opzuigpositie kan de positie storen en kan dan leiden tot onverwachte fouten (bijv. **DiTi verloren**).

 Zorg ervoor dat er geen sterke magneet in een SBS-positie naast de aanzuigpositie wordt geplaatst.



Fluent ID-segmenten kunnen alleen worden verwijderd door de technicus vanwege de aansluiting op de elektroplaat onder het dek.

- ✓ Methode moet door de key-operator worden voorbereid.
- De verbruiksmiddelen komen overeen met de verbruiksmiddelen die in de methode zijn gedefinieerd.
- Fluent ID-buisrunners mogen alleen worden geladen nadat de methode werd gestart, als dit wordt gevraagd op het touchscreen.
- 1. Volg de op het touchscreen weergegeven instructies op.





De afbeelding toont een voorbeeld van een instructie (A) die op het touchscreen wordt weergegeven:

### 6.4 Alvorens een methode te starten

De volgende checklist moet worden voltooid voordat een methode wordt gestart.

Instrument/component	Taak	Referenties/activiteiten
Procesvalidatie	Controleer voor het starten van een produc- tiecyclus of de geselec- teerde methode werd gevalideerd.	Raadpleeg de key-ope- rator voor meer infor- matie.
Touchscreen	Volg de instructies op het touchscreen op. LET OP! Deze instructies van de key-operator moeten strikt in acht worden genomen.	_
	Als er geen instructies verschijnen, de onder- staande takenlijst op- volgen.	

Tab. 15: Controles voor het starten van een methode



Instrument/component	Taak	Referenties/activiteiten
Segmenten, dragers, opties en apparaten	Zorg ervoor dat alle segmenten, dragers, opties en apparaten zijn geïnstalleerd en bevei- ligd. Zorg ervoor dat alleen objecten die bij deze methode worden ge- bruikt op het dek aan- wezig zijn. Zorg ervoor dat de test- cyclus succesvol wordt afgesloten.	Als de testcyclus mis- lukt, contact opnemen met de key-operator om de testcyclus op- nieuw uit te voeren.
Monsters en reagentia	Zorg ervoor dat alle monsters, reagentia en al het laboratoriumma- teriaal correct zijn gela- den. LET OP! Het scannen van barcodes wordt pas uitgevoerd nadat de methode is gestart. Zorg ervoor dat er zich op het Fluent ID-dek geen runners bevinden, alvorens de methode te starten. De runners mo- gen alleen worden gela- den, als dit wordt ge- vraagd op het touchs- creen.	
Afvoerbuizen (alleen vloeistofsystemen)	Zorg ervoor dat de af- voerbuis correct ver- loopt.	Voer een visuele con- trole uit van de afvoer- buizen om er zeker van te zijn dat deze niet ge- knikt of geplet zijn. Vervang defecte afvoer- buizen. Zie paragraaf "Wasstation aansluiten (MCA 384)" [▶ 170].
Wassysteem (alleen vloeistofsystemen)	Zorg ervoor dat sys- teemvloeistof en af- voerreservoir correct zijn verbonden.	Zie paragraaf "Buizen van systeemvloeistofre- servoir en afvoerreservoir controleren" [> 93].



Instrument/component	Taak	Referenties/activiteiten		
Wassysteem (alleen vloeistofsystemen)	Zorg ervoor dat het sys- teemvloeistofreservoir tot de correcte stand is gevuld. Zorg ervoor dat het af- voerreservoir leeg is.	Zie paragraaf "Het sys- teemvloeistofreservoir en afvoerreservoir aanslui- ten" [▶ 146].		
Wassysteem (alleen vloeistofsystemen)	Zorg ervoor dat de cor- recte systeemvloeistof zoals gedefinieerd in de methode wordt ge- bruikt.	_		
Wassysteem (alleen MCA-wassysteem)	Controleer het vloeistof- niveau in het wasblok.	-		
Wegwerppunt-afvoer en wasstationeenheid	Zorg ervoor dat de wegwerppunt-afvoer en de wasstationeenheid schoon zijn. Zie paragraaf "We werppunt-afvoer en stationeenheid rein [▶ 140].			
	Zorg ervoor dat afdek- kingen voor afvoergoten voor aerosol-inperking of voor MCA 384-punt- geleiding zijn gemon- teerd.	_		
Wegwerppunten	Zorg ervoor dat de cor- recte punten worden geladen. Zorg ervoor dat de pun- tafvoer leeg is.	-		
Vaste punten	Zorg ervoor dat de vas- te punten schoon en onbeschadigd zijn.	Voer een visuele con- trole op reinheid van de vaste punten uit. Voer met een tandarts- spiegel een visuele controle op correcte		
		toestand van de coating uit.		
Dek	Zorg ervoor dat op het dek geïnstalleerde dra- gers, laboratoriumma- teriaal en apparaten overeenkomen met de dekindeling die voor deze methode is gede- finieerd.	Zie paragraaf "Dekinde- ling controleren" [▶ 89].		



Instrument/component	Taak	Referenties/activiteiten
Laboratoriummateriaal	Zorg ervoor dat al het laboratoriummateriaal correct is gepositio- neerd. Als de multititer- platen zijwaartse bewe- ging vertonen, moet er- voor worden gezorgd dat de positioneervoor- zieningen voor labora- toriummateriaal correct zijn.	Zie paragraaf "Laborato- riummateriaal positione- ren" [▶ 171].
Buisrotator	Controleer of er geen positioneerpennen of vergrendelpennen op de buisrotator ontbreken.	Zie paragraaf "Vergren- delpennen en positio- neerpennen vervangen" [▶ 198]

### 6.4.1 Buizen van systeemvloeistofreservoir en afvoerreservoir controleren

### **▲ VOORZICHTIG**

### Monsterverontreiniging!

Bij een dubbele vloeistof-configuratie van de FCA kunnen verschillende systeemvloeistoffen voor elke arm worden gebruikt. Aansluiten van het onjuiste systeemvloeistofreservoir op een arm kan leiden tot verontreiniging van monsters.

- Label elk systeemvloeistofreservoir met de bijbehorende systeemvloeistofnaam.
- ✓ Het wassysteem moet correct worden geïnstalleerd.
- 1. Controleer of buis (B) correct met het systeemvloeistofreservoir (D) is verbonden.





2. Indien aanwezig, controleren of vloeiststofdetectiesensor (C) correct is verbonden met het systeemvloeistofreservoir (D).

### 6.4.2 Buizen van afvoerreservoir controleren

1. Controleer of buis (B) correct met het afvoerreservoir (D) is verbonden.



- 2. Controleer of de buizen (C) en (D) correct zijn verbonden met het afvoerreservoir.
- 3. Indien aanwezig, controleren of vloeiststofdetectiesensor (E) correct is verbonden met het afvoerreservoir.



4. Schroef deksel (F) erop.



### 6.5 Een methode uitvoeren

Een methode is een verzameling scripts of processen zoals gedefinieerd in de FluentControl-software. De methode kan worden uitgevoerd in een cyclus.

De key-operator schrijft een methode die als volgt kan worden uitgevoerd.

### LET OP

### Instrumentbeschadiging!

As het dek niet correct is ingesteld of als de software niet correct werkt of verkeerd wordt gebruikt, kan dit leiden tot instrumentbeschadiging.

- Zorg ervoor dat alle veiligheidsvoorzieningen zijn geïnstalleerd en functioneren.
- Zorg ervoor dat op het dek geïnstalleerde dragers, laboratoriummateriaal en apparaten overeenkomen met de dekindeling die voor deze methode is gedefinieerd.
- Zorg ervoor dat alleen objecten die bij deze methode worden gebruikt op het dek aanwezig zijn.

### 6.5.1 Een methode starten

- ✓ FluentControl is gestart.
- ✓ Paragraaf "Gebruikersaanmelding" [▶ 83] is uitgevoerd.
- ✓ Paragraaf "Alvorens een methode te starten" [▶ 90] is uitgevoerd.
- 1. Selecteer Method Starter (Methodestarter) (A).





Toets wordt verlicht zodra deze wordt aangeraakt.

- 2. Selecteer de methode (B) die moet worden uitgevoerd. *Geselecteerde methode is gemarkeerd.*
- 3. Druk op OK.





4. Druk op Run (Uitvoeren) (C).



- 5. Volg de instructies op het touchscreen op.
- Als uw script DeckCheck omvat, let dan op verschillen van de momentele livedekindeling vergeleken met de verwachte referentie-dekindeling. Zie "DeckCheck-bedrijf" [▶ 107].
- 7. Wacht tot de methodecyclus is beëindigd.

Het scherm toont de geschatte tijd (D) waarop de methode zal eindigen.







8. Als het bericht **Run finished with error(s)!** (Run uitgevoerd met fout(en)!) wordt weergegeven, dan drukt u op **Log** (Logboek) om de fouten en waarschuwingen te controleren.

TouchTools	Running 🛞	0
1		
	Run finished with error(s)	!
	• •	
	11:59 Start Time	12:19 End Time

9. Druk op Next (Volgende) om terug te keren naar het beginscherm.

Ru	n finished v	vith er	ror(s)!
11:59 Start Time			12: End Ti
Comment	Process	Log	
2 E: Script exec 1 Run paused.	ution terminated by End Script st.	atement.	12:00:48 (2017-08-18) 11:59:35 (2017-08-18)

10. Druk op **Exit** (Afsluiten).





Het scherm hieronder (E) wordt weergegeven, als de methodecyclus eindigt.

6.5.2 Fluent ID-runners laden en ontladen



### De Fluent een klasse 1 laserproduct conform IEC 60825-1:2014 dat laserstraling uitzendt.

Verblinding, flitsblindheid en nabeelden kunnen door de laserstraal worden veroorzaakt.

Staar niet in de laserstraal of in de reflecties ervan.

#### 6.5.2.1 Fluent ID-runners laden

### LET OP

#### Beschadiging door verkeerd in- en uitladen

Schade aan runners en pennen.

- Lijn de runner horizontaal uit ten opzichte van het dek.
- Ondersteun het voorste einde van de runner met een hand.
- Zorg er tijdens uitladen voor dat de runner alle pennen vrijmaakt, alvorens de runner op te tillen.
- ✓ Fluent is uitgerust met een Fluent ID-buisbarcodescanner.
- ✓ Buizen worden geladen in de runners met het barcodelabel naar links.
- ✓ Alle buizen in een runner hebben dezelfde afmeting en vorm. Voor buisrunnertypen, zie paragraaf "Met Fluent ID compatibele buisrunners" [▶ 61].
- 1. Selecteer en start de methode met het touchscreen.

De leds knipperen en de melding **Buizen laden** wordt weergegeven op het touchscreen.



Bij gebruik van verschillende soorten buizen moet ervoor worden gezorgd dat voor elk specifiek raster het juiste type runner wordt gebruikt.

- 2. Ondersteun het voorste einde van de runner met een hand.
- 3. Houd de runner horizontaal ter hoogte van het dek.
- 4. Duw de runner tegen de stoppositie.



5. Schuif de buisrunners achtereenvolgens op de rasters in het Fluent IDlaadbereik.

₽	Running	Load tube r	unn	ers								
			000000000000000000000000000000000000000		000000000000000000000000000000000000000	4	000000000000000000000000000000000000000	6				
	Ple	ase load tube runners and	pres	s cont	tinue	when	finish	ed		Errors:	] Missin	a Runner
										쎚		~
										Stop		Continue

6. Controleer of alle barcodes succesvol werden gescand. De leds branden groen, als de buisrunners in geladen positie zijn en alle barcodelabels succesvol werden gescand.



Voor de beschrijving van Fluent ID-ledstatus, zie paragraaf "Fluent IDstatusleds" [> 79].



- 7. In geval van een barcodescanfout, de runner ontlasten, het probleem oplossen en de runner weer beladen.
- 8. Trek de runner horizontaal tot volledige verwijdering langs het dek.



De Fluent ID leest elke code meerdere keren, als deze de scanner passeert. Voor kleine en smalle buizen (bijv. diameter  $\leq$  10 mm) de snelheid van handmatig laden reduceren om alle lees- en foutmeldingen te beperken.

00000000000		000000000000000000000000000000000000000	0000000000000	000000000000000000000000000000000000000	000000000000000000000000000000000000000											
1	2	3	4	5	6											
Errors:		Missi	ng Ru	nner	D V	Vrong	g Run	ner	•	Unrea	dable	•	Duplicate	s (	С	Missing

Errors. - Missing Kumer - Missing Kumer - Sineddable - Supileates - Missin

Afb. 42: Barcodeleesbevestiging weergegeven op het touchscreen

#### Tab. 16: GUI betekenis (runner)

Vierkant (runner)	Betekenis
Groen	Alle buisbarcodes in runner succesvol gelezen.
Wit met rode contour	Onjuist type runner voor deze raster- positie.



Vierkant (runner)	Betekenis
Grijs met rode contour	Ontbrekende runner. Een runner moet in deze rasterpositie worden geladen.

#### Tab. 17: GUI betekenis (buispositie)

Cirkel (buispositie)	Betekenis
Groen	Barcodes succesvol gelezen.
Rood	Onleesbare barcode
Oranje	Barcode dupliceren
Wit met rode contour	Ontbrekende buis. Een buis moet in deze positie worden geladen.

6

Als de 2 ml Safe-Lock-buisrunner wordt gebruikt, is het niet mogelijk om ontbrekende buizen te onderscheiden van onleesbare barcodes. Ontbrekende buizen worden gerapporteerd als niet-leesbare barcodes.

#### 6.5.2.2 Fluent ID-runners uitladen

- ✓ De cyclus is voltooid of er is een cyclus actief en de leds knipperen met de melding Buizen ontladen weergegeven op het touchscreen.
- 1. Trek de runner horizontaal tot volledige verwijdering langs het dek.

#### 6.5.3 Buisrotator-runners in- en uitladen

#### 6.5.3.1 Buisrotator-runners inladen

### **A VOORZICHTIG**

#### Biologische verontreiniging van het systeem en/of gebruiker!

Beschadigde monsterbuizen kunnen imploderen; dit kan leiden tot morsen van het monster op de buisrotator.

- · Voorkom dat beschadigde buizen op de buisrotator worden geladen.
- ✓ De Fluent is uitgerust met een buisrotator.
- Buizen worden ingeladen in de buisrotator-runners met het barcodelabel naar links.
- ✓ Alle buizen in een runner hebben dezelfde afmeting en vorm. Voor buisrunnertypen, zie paragraaf "Buisrotator-runners" [▶ 64].
- 1. Selecteer en start de methode met het touchscreen.

De leds knipperen en de melding **Buizen laden** wordt op het touchscreen weergegeven.

Bij gebruik van verschillende buistypen, ervoor zorgen dat het correcte runnertype voor het buistype wordt gekozen (BD of Greiner). Zorg ervoor dat



verschillend hoge buizen op de bijbehorende runners worden geladen: De buizen worden altijd door de runnerbrug ter hoogte van hun doppen op hun plaats gehouden. De buisonderzijde moet altijd correct in de buisinzetstukken van de runners worden geplaatst.

2. Open de runner-vergrendelingshendel.



3. Ondersteun het voorste einde van de runner met een hand.



- 4. Houd de runner horizontaal ter hoogte van het dek.
- 5. Duw de runner tegen de stoppositie.



$m{\Delta}$ $ angle$ Running $ angle$ Load tube runners	
Please load tube runners and press continue when finished	Errors: 🔲 Missing Runner

6. Schuif de buisrotator-runners achtereenvolgens op de rasters van de buisrotator.

7. Controleer of alle barcodes succesvol werden gescand.

De leds branden groen, als de buisrunners in geladen positie zijn en alle barcodelabels succesvol werden gescand. Voor de beschrijving van de buisrunner-ledstatus, zie paragraaf "Fluent IDstatusleds" [> 79].

- 8. In geval van een barcodescanfout, de runner ontlasten, het probleem oplossen en de runner weer beladen.
- 9. Sluit de runner-vergrendelingshendel.







Afb. 43: Barcodeleesbevestiging weergegeven op het touchscreen

#### Tab. 18: GUI betekenis (runner)

Vierkant (runner)	Betekenis
Groen	Alle buisbarcodes in runner succesvol gelezen.
Wit met rode contour	Onjuist type runner voor deze raster- positie.
Grijs met rode contour	Ontbrekende runner. Een runner moet in deze rasterpositie worden geladen.

#### Tab. 19: GUI betekenis (buispositie)

Cirkel (buispositie)	Betekenis
Groen	Barcodes succesvol gelezen.
Rood	Onleesbare barcode
Oranje	Barcode dupliceren
Wit met rode contour	Ontbrekende buis. Een buis moet in deze positie worden geladen.

### 6.5.3.2 Buisrotator-runner uitladen

6

Bewaar geen buisrotator-runners die ingeladen zijn met buizen buiten de bedrijfsomstandigheden van de vloeistofverwerking. Zie paragraaf "Omgevingsvoorwaarden" [> 45].

- De cyclus is voltooid of er is een cyclus actief en de leds knipperen met de melding Buizen ontladen weergegeven op het touchscreen.
- ✓ De buisrotator is in horizontale uitgangspositie.



1. Open de runner-vergrendelingshendel.



2. Trek de runner horizontaal tot volledige verwijdering langs het dek.



Buisrotators kunnen alleen door de technicus worden verwijderd vanwege de aansluiting op de elektroplaat onder het dek.

- Methode moet door de key-operator worden voorbereid.
- De verbruiksmiddelen komen overeen met de verbruiksmiddelen die in de methode zijn gedefinieerd.
- Buisrotator-runners mogen alleen worden ingeladen nadat de methode werd gestart, als dit wordt gevraagd op het touchscreen.

#### 6.5.4 Fouten resetten

Als een melding wordt weergegeven, als volgt te werk gaan:

- Melding
- 1. Controleer de werking van het display, de toets of foutmeldingen. Zie paragrafen "Werkbereik" [▶ 73] en "Methodeherstel-toetsen" [▶ 74].
- 2. Volg de instructies in deze handleiding en op het touchscreen om de fout te corrigeren.
- 3. Ga verder met de methodecyclus. Zie paragraaf "Display, optie en actietoetsen" [▶ 73].

Als de statuslamp gaat branden of als deze van kleur verandert, als volgt te werk gaan:

Statuslamp

- 1. Controleer de instrumentstatus. Zie paragraaf.
- Als de Fluent is uitgerust met een Fluent ID-buisbarcodescanner, de ledstatus van de Fluent ID-buisbarcodescanner controleren. Zie paragraaf "Fluent IDstatusleds" [▶ 79].
- 3. Controleer de werking van het display, de toets of foutmeldingen. Zie paragrafen "Werkbereik" [▶ 73] en "Methodeherstel-toetsen" [▶ 74].



- 4. Controleer de tabel voor het verhelpen van storingen. Zie paragraaf "Tabellen voor het verhelpen van storingen" [▶ 157].
- 5. Als het probleem niet kan worden opgelost, contact opnemen met "Klantsupport" [▶ 203].

### 6.6 DeckCheck-bedrijf

Als uw script gebruik van de DeckCheck omvat, maakt het DeckCheckcamerasysteem afbeeldingen van de werktafel na het laden om deze liveafbeeldingen met de referentie te vergelijken.

De DeckCheck heeft ongeveer 20 seconden nodig voor een 3-armig systeem / 3 camera's en ongeveer 12 seconden voor een systeem met 1 of 2 armen/ enkele camera om afbeeldingen van het dek te maken en de vergelijking tussen de liveen referentie-dekindelingen weer te geven (ervan uitgaande dat de pc-configuratie geschikt is – zie de handleiding voor toepassingssoftware FluentControl).

Bij de eerste keer dat de DeckCheck-opdracht wordt gebruikt nadat het instrument is ingeschakeld, kan het enkele minuten duren voordat het eerste resultaat wordt weergegeven.

Gedurende deze tijd wordt de achter-LED ingeschakeld.

Bij 3-armige Fluent-systemen moet de middenarm tussen de linker- en rechterpositie bewegen (bij 1- of 2-armige systemen zullen de linker- en rechterarm respectievelijk uiterst links en uiterst rechts worden geplaatst.) Voor deze armbeweging moet de voordeur gesloten zijn. Als de afbeelding wordt gemaakt terwijl de deur open is, wordt één camera geblokkeerd door de middenarm of door een arm die handmatig werd bewogen.

Tijdens het DeckCheck-proces beeldt het touchscreen vormen af die over het scherm bewegen en **Afbeeldingen maken** gevolgd door **Controleren**. Na 12–20 seconden, afhankelijk van de instrumentgrootte en de configuratie worden dekafbeeldingen afwisselend weergegeven. De weergegeven afbeelding zal de eerste camera zijn met een geconstateerde afwijking, beginnend vanaf links.

DeckCheck-scherm geeft een afwijking ten opzichte van de referentie-afbeelding. Hier toont de referentie-afbeelding dat er een plaat aanwezig zou moeten zijn en dat de middencamera de afwijking heeft gedetecteerd.



Als de **Ignore & Continue** toets verschijnt, wordt de cyclus voortgezet met de actuele werktafel. Selecteer de **Ignore & Continue** toets, als u er zeker bent dat er geen verschillen meer zijn met de vereiste werktafel, en sluit dan de deur. Select



**Check** als u nieuwe afbeeldingen van het dek wilt maken — neem in acht dat, indien de deur bij een 3-armig systeem niet gesloten is, de afbeelding wel wordt opgenomen maar dat de middenarm de camera zal blokkeren. Anders wordt automatisch een controle uitgevoerd, als de deur gesloten is.



Als het script de optie **altijd weergeven**bevat, wordt bovenstaand scherm weergegeven en worden er geen afwijkingen weergegeven. De referentieafbeeldingen en live-afbeeldingen wisselen elkaar echter af en het kan zijn dat er kleine afwijkingen zijn die niet door het systeem worden geregistreerd, maar die met het menselijke oog eenvoudig kunnen worden waargenomen – bijvoorbeeld kleurverschillen, enkele ontbrekende buizen/tips of kleine zijdelingse verschuivingen. Zie de hieronder vermelde grenswaarden.

Als afwijkingen worden waargenomen, worden deze gemarkeerd.

Afwijkingen corrigeren:

- 1. Open de deur en vervang of corrigeer de positie van de gemarkeerde delen.
- 2. De DeckCheck vergelijkt de gecorrigeerde live-situatie ononderbroken met de referentie-dekindeling.
- Bekijk met de DeckCheck-toetsen de afwijkingen zoals waargenomen door elk van de camera's of pauzeer de afbeelding en vergelijk deze indien nodig met de referentie-afbeelding. *Als er geen afwijkingen meer worden waargenomen, verschijnt de groene Continue-toets.*
- 4. Select **Continue** om de methode voort te zetten.






Als resterende verschillen als acceptabel worden beoordeeld (bijvoorbeeld het totaal aantal tips varieert aan het begin van de methode, of vloeistofniveaus wijken aanzienlijk af aan het begin van een cyclus), kunt u **Ignore & Continue** selecteren, als dit in het script van uw key-operator wordt weergegeven.

Sommige afwijkende dekindelingen kunnen niet worden gemarkeerd door de DeckCheck—bijvoorbeeld de volgende gekleurde FCA-tip-trays:

Verschil tussen typen MCA-kopadapter:

- Geel/oranje
- Wit/oranje
- Grijs/alle kleuren

MCA 384 verschillende tip-typen

MCA 96 verschillende tip-typen

Ontbrekende buizen op gedeeltelijk geladen buisrunners

Bak 300 SBS

Microplaten 180 graden gedraaid

Microtiterplaat well-vormig (bijvoorbeeld ronde versus platte bodem of PCR-well)

Platen in perifere hotels 10 ml/ 25 ml bakken als inzetstuk

Sommige transparante deksels

Sommige van deze verschillen zijn echter duidelijk zichtbaar bij de schakeling tussen live- en referentie-dekindelingen.

### 6.7 Methodeherstel

FluentControl biedt de optie om van fouten te herstellen-bijv .:

Eerdere methode werd afgebroken of er is een fatale fout opgetreden: De optie voor methodeherstel biedt de mogelijkheid tot voortzetten vanaf het punt waarop deze fout tijdens een eerdere cyclus is opgetreden.



Nadat een methode werd afgebroken of er een fatale fout is opgetreden, moet dagelijks onderhoud worden uitgevoerd. Zie paragraaf "Dagelijkse systeemverzorging" [> 117].

#### 6.7.1 Wisselen naar methodeherstelmodus

- ✓ Key-operator heeft de optie voor methodeherstel in FluentControl geactiveerd.
- ✓ Eerdere methodecyclus werd afgebroken.



1. Selecteren Methodeherstel (A).



### 6.7.2 Een methodecyclus herstellen

- ✓ Paragraaf "Wisselen naar methodeherstelmodus" [▶ 109] werd uitgevoerd.
- 1. Selecteer verder naar het volgende scherm (A).

Het scherm toont de laatst uitgevoerde scriptregel (C) en de scriptregel waarbij de fout is opgetreden—het herstelpunt (B).



- 2. Selecteer vereiste toetsen die zijn beschreven in paragraaf "Methodehersteltoetsen" [▶ 74].
- 3. Zorg ervoor dat de fysische dekindeling van de Fluent overeenkomt met de werktafelindeling (D) die wordt weergegeven op het touchscreen.
- 4. Selecteren Cyclusherstel (E).



### Het systeem start.



# 6.8 Het instrument uitschakelen

Als er geen methode actief is, wisselt het instrument naar stand-bymodus. Het instrument moet niet van het net worden losgekoppeld.

Schakel het instrument als volgt uit:

- 1. Stop alle methodes en selecteer de stand-bymodus op het touchscreen.
- 2. Zorg ervoor dat het instrument in stand-bymodus is. Zie paragraaf.
- 3. Plaats de robotarmen in een bewegingsvrij bereik.





4. Schakel de voeding uit met voedingsschakelaar (A) aan de achterzijde van de externe stroomvoorziening.



# 7 Systeemverzorging

Dit hoofdstuk geeft instructies over alle taken die moeten worden uitgevoerd om de Fluent in goede toestand te houden.



Gebruik de Fluent uitsluitend, als deze in goede werktoestand is. Neem instructies voor systeemverzorging zoals in deze handleiding beschreven in acht. Om optimale prestaties en betrouwbaarheid te garanderen, de onderhouds- en reinigingstaken regelmatig uitvoeren.

Raadpleeg bij eventuele problemen en vragen de paragraaf "Klantsupport" [> 203].



# A VOORZICHTIG

#### Arm botst tegen werkbare objecten aan

Vloeiende armen kunnen handmatig worden bewogen. Zorg ervoor dat de handmatige bewegingen van de armen soepel worden uitgevoerd terwijl de arm wordt vastgehouden. Stoot de armen niet tegen vaste voorwerpen, waaronder de mechanische stop voor het bereik van de armen

# 7.1 Ontsmetting

Ontsmetting is volgens de standaard laboratoriumregelgeving vereist onder de omstandigheden zoals beschreven in paragraaf "Ontsmettingsverklaring" [> 39].

### **A WAARSCHUWING**

### Verontreiniging!

Resten van stoffen op de Fluent kunnen leiden tot letsel en kunnen de integriteit van het proces nadelig beïnvloeden.

 Ontsmet de Fluent en zijn onderdelen en accessoires vóór alle andere werkzaamheden.

De ontsmettingsmethode moet zijn gedefinieerd door de key-operator, gebaseerd op het type en de mate van verontreiniging. Richtlijn bij de selectie van de ontsmettingsmiddelen en toepassingsmethodes staan in dit hoofdstuk.



Voor informatie over de behandeling met waterstofperoxide-dampen, zie de referentiehandleiding. Zie Referentiedocumenten.

# 

### Onjuiste meetresultaten van de Frida Reader!

Zonder gemonteerd inzetstuk kan de Frida Reader onjuiste meetresultaten leveren.

 Gebruik de rode, blinde stop, als het inzetstuk wordt verwijderd (bijv. voor reiniging).



# 7.2 Reinigingsmiddelen

### 7.2.1 Reinigingsmiddelen specificaties

Speciale reinigingsmiddelen zijn voor de systeemverzorging vereist. Alle aanbevolen reinigingsmiddelen werden zorgvuldig geselecteerd en getest.

### LET OP

### Gereduceerde effectiviteit en chemische compatibiliteit!

Er is geen garantie voor de effectiviteit van reinigingsmiddelen en de chemische compatibiliteit, als andere reinigingsmiddelen dan die aanbevolen door Tecan worden gebruikt.

- Gebruik alleen reinigingsmiddelen die door Tecan worden aanbevolen.
- Reinigingsmiddelen worden in de tabellen voor systeemverzorging voor elke specifiek gebruik gedefinieerd. Gebruik geen reinigingsmiddelen, als deze niet voor een speciale taak zijn gespecificeerd.

In de volgende tabel staan de reinigingsmiddelen die zijn gespecificeerd voor gebruik zoals beschreven in de onderhoudstabellen van het systeem en de onderhoudsactiviteiten van het systeem:

*Tab. 20:* Reinigingsmiddelen voor gebruik zoals beschreven in de onderhoudstabellen van het systeem

Middel	Specificatie
DI-water	Gedestilleerd of gede-ioniseerd wa- ter
Alcohol	70% ethanol, 100% isopropanol (2-propanol)
Mild reinigingsmiddel	Liqui-Nox
Oppervlaktereinigingsmiddel	Contrad 70, Contrad 90 / Contrad 2000, Decon 90
Ontsmettingsmiddel	Bacillol plus, SporGon
Ontsmettingsmiddel voor oppervlak- ken (voor besmetting met nucleïnezuur)	DNAzap
Mild zuur	zwavelzuur 0,3M, 10% azijnzuur, 30-40% mierenzuur
Base	Natriumhydroxide 0.1M
Bleekmiddel	2% natriumhypochloriet



Middel	Specificatie
Systeemvloeistof	Zoals gedefinieerd in de methode. Neem in acht dat waterige oplossin- gen met zoutgehalte moeten worden uitgespoeld, als het systeem niet ac- tief is-bijvoorbeeld 's nachts of in het weekend. Zie Systeemverzorging "Einde van de dag" [> 119].

### 7.2.2 In de handel verkrijgbare reinigingsmiddelen

Alle instructies-door de fabrikant van de reinigingsmiddelen gegeven of in deze handleiding-voor de omgang met reinigingsmiddelen moeten zorgvuldig worden gelezen en opgevolgd.

In de onderstaande tabel staan een aantal in de handel verkrijgbare reinigings- en desinfectiemiddelen, gespecificeerd voor gebruik zoals beschreven in de onderhoudstabellen van het systeem en de onderhoudsactiviteiten van het systeem.

Reinigingsmiddel	Middelcategorie	Fabrikant
DNAzap	Ontsmettingsmiddel voor oppervlakken (voor oppervlakken verontrei- nigd met nucleïnezuren)	Ambion www.ambion.com
Decon, Contrad	Oppervlaktereinigings- middel	Decon Laboratories www.deconlabs.com
SporGon	Ontsmettingsmiddel	Decon Laboratories www.deconlabs.com
Bacillol Plus	Ontsmettingsmiddel	www.bode-chemie.com
Liqui-Nox	Mild reinigingsmiddel	Alconox www.alconox.com

Tab. 21: In de handel verkrijgbare reinigingsmiddelen

## 7.3 Systeemverzorgingsmodus

De key-operator definieert de vereiste systeemverzorgingsmethoden zoals weergegeven in systeemverzorgingstabellen in "Systeemverzorgingstabellen" [▶ 117]. Het **Systeemverzorging** op het touchscreen biedt oriëntatie voor systeemverzorgingstaken.

### 7.3.1 Naar systeemverzorgingsmodus wisselen

✓ Er moeten systeemverzorgingsmethodes beschikbaar zijn.



- A System Care Method Starter E
- 1. Selecteren **Systeemverzorging** (A).

2. Selecteer de taak die moet worden uitgevoerd.



- 3. Druk op Cyclus om de systeemverzorgingsmodus te initiëren.
- 4. Voer de systeemverzorgingstaken uit.

### 7.3.2 Fouten resetten

Als een melding wordt weergegeven, als volgt te werk gaan:

Melding

- 1. Controleer de werking van het display, de toets of foutmeldingen. Zie paragrafen "Werkbereik" [▶ 73] en "Methodeherstel-toetsen" [▶ 74].
- 2. Volg de instructies in deze handleiding en op het touchscreen om de fout te corrigeren.
- 3. Ga verder met de methodecyclus. Zie paragraaf "Display, optie en actietoetsen" [▶ 73].



Als de statuslamp gaat branden of als deze van kleur verandert, als volgt te werk gaan:

#### Statuslamp

- 1. Controleer de instrumentstatus. Zie paragraaf.
- Als de Fluent is uitgerust met een Fluent ID-buisbarcodescanner, de ledstatus van de Fluent ID-buisbarcodescanner controleren. Zie paragraaf "Fluent IDstatusleds" [▶ 79].
- 3. Controleer de werking van het display, de toets of foutmeldingen. Zie paragrafen "Werkbereik" [▶ 73] en "Methodeherstel-toetsen" [▶ 74].
- 4. Controleer de tabel voor het verhelpen van storingen. Zie paragraaf "Tabellen voor het verhelpen van storingen" [▶ 157].
- 5. Als het probleem niet kan worden opgelost, contact opnemen met "Klantsupport" [▶ 203].

### 7.4 Systeemverzorgingstabellen

Om optimale prestaties en betrouwbaarheid te garanderen, de onderhouds- en reinigingstaken uitvoeren zoals aanbevolen.



De taken in de systeemverzorgingstabel kunnen alleen worden uitgevoerd in de systeemverzorgingsmodus. Zie paragraaf "Systeemverzorgingsmodus" [> 115].

De systeemverzorgingstaken moeten worden uitgevoerd in regelmatige intervallen —namelijk dagelijkse, wekelijkse en maandelijkse systeemverzorging.

#### 7.4.1 Dagelijkse systeemverzorging

#### 7.4.1.1 Begin van de dag

Voer de methode **DailySystemCare** methode uit, indien beschikbaar gesteld door de key-operator, of voer elke individuele taak uit die van toepassing is op uw Fluent-armconfiguratie, in chronologische volgorde vermeld in de onderstaande tabel.

Tab. 22: Systeemverzorgingstabel voor begin van de dag

Instrument/ component	Systeemverzorgings- taak	Reinigingsmiddel/ wegwerpproduct/ apparaat	Referentie/ systeemverzorgingsactivi- teiten
Doorprikpunten	Controleer de door- prikpunten visueel op eventuele afzettingen. Indien nodig reinigen. Controleer of de pun- ten verbogen zijn.	70% ethanol of 2% bleekmiddel en pluis- vrije doek	Zie paragraaf <b>"Doorprikpun-</b> ten reinigen" [▶ 131].



Instrument/ component	Systeemverzorgings- taak	Reinigingsmiddel/ wegwerpproduct/ apparaat	Referentie/ systeemverzorgingsactivi- teiten
Wegwerppunt-kegels en vaste naalden	Op beschadiging en afzettingen controle- ren	_	Deze taak maakt deel uit van de <b>dagelijkse systeem-</b> verzorgings uit.
			LET OP! Wegwerpnaalden zijn niet bedoeld voor her- gebruik.
Vaste naalden	Reinigen. Controleer of de punten verbo- gen zijn. Controleer visueel met een tand- artsspiegel of de coating intact is.	70% ethanol of 100% isopropanol en pluis- vrije doek	Zie paragraaf "Vaste punten reinigen" [▶ 131].
Systeemvloeistofre- servoir (Liquid FCA en MCA met vaste paalden)	Ervoor zorgen dat de- ze schoon en vol is, zonder zichtbare bel- len is	_	Deze taak maakt deel uit van de <b>dagelijkse systeem-</b> verzorgings uit.
	Ervoor zorgen dat de slangen naar de re- servoirconnectors correct zijn aangeslo- ten		
Vloeistof-afvoerreser- voir (Liquid FCA en MCA met vaste naalden)	Ervoor zorgen dat de- ze leeg is Ervoor zorgen dat de slangen naar de re- servoirconnectors correct opnieuw zijn aangesloten	-	Deze taak maakt deel uit van de <b>dagelijkse systeem-</b> verzorgings uit.
Wegwerppunt-afvoer- zak	Ervoor zorgen dat de- ze leeg is	_	Zie paragraaf "Wegwerp- punt-afvoerzak vervangen" [▶ 143].
			Deze taak maakt deel uit van de <b>dagelijkse systeem-</b> verzorgings uit.
Vloeistofsysteem (Liquid FCA)	Ervoor zorgen dat de- ze schoon is	Systeemvloeistof, al- cohol, DI-water	Deze taak maakt deel uit van de <b>dagelijkse systeem-</b> verzorgings methode of kan apart worden uitgevoerd als Liquid FCA routinematig spoelonderhoud uit. Zie paragraaf "Vloeistofpad
			reinigen" [> 145].



Instrument/ component	Systeemverzorgings- taak	Reinigingsmiddel/ wegwerpproduct/ apparaat	Referentie/ systeemverzorgingsactivi- teiten
Wassysteem (MCA met vaste naal- den)	Spoelen/primen	Wasvloeistof	De directe opdracht uitvoe- ren <b>Wasstation primen</b> (MCA384)uit.
Vloeistofsysteem (Liquid FCA)	Visueel controleren op afwezigheid van drup- pels op de punten of DiTi-kegel na spoelen	-	Deze taak maakt deel uit van de <b>dagelijkse systeem-</b> verzorgings uit.
Grijpervingers	Controleren of vingers recht en waterpas zijn Controleren op be- schadiging en onjuiste uitlijning	_	Bij onjuiste uitlijning, zie pa- ragraaf "Robotgrijperarm (RGA) verhelpen van storin- gen" [▶ 166]. Vervorming of beschadi- ging. Raadpleeg de para- graaf "Klantsupport" [▶ 203].
FCA-grijpervingers	Op beschadiging controleren	-	Vervangen indien bescha- digd. Voor de bestelgege- vens, zie de referentiehand- leiding. Zie de paragraaf Referentiedocumenten.
Frida Reader	Verwijder de blinde plug en steek het in- zetstuk in de Frida Reader	-	Zie paragraaf "Frida Reader" [▶ 153].

# **⚠ VOORZICHTIG**

### Onjuiste meetresultaten van de Frida Reader!

Zonder gemonteerd inzetstuk kan de Frida Reader onjuiste meetresultaten leveren.

• Gebruik de rode, blinde stop, als het inzetstuk wordt verwijderd (bijv. voor reiniging).

### 7.4.1.2 Einde van de dag

De volgende tabel somt in chronologische volgorde de dagelijkse systeemverzorgingstaken aan het einde van de dag op:



### Tab. 23: Systeemverzorgingstabel voor eind van de dag

Instrument/	Systeemverzorgings-	Reinigingsmiddel/	Referentie/	
component	taak	wegwerpproduct/ apparaat	systeemverzorgingsactivi- teiten	
Dektrays	Controleer op mor- sen. Reinigen of ver- vangen indien nodig.	DI-water, alcohol, mild reinigingsmiddel, ontsmettingsmiddel, base, bleekmiddel, DNAzap	Zie paragraaf "Dektrays rei- nigen" [▶ 136].	
Segmenten Fluent ID-behuizing	Reinigen	DI-water, alcohol, mild reinigingsmiddel, ontsmettingsmiddel, base, bleekmiddel, DNAzap	Zie paragraaf "Runners and segmenten reinigen" [> 136]. WAARSCHUWING! Kijk niet in de laserstraal.	
		venster vereist andere reinigingsmiddelen dan het segment zelf. Zie "Wekelijkse sys- teemverzorging" [▶ 123].		
Reflectorfolie (Fluent ID, buisrota- tor)	Reinigen en op be- schadiging controle- ren	Alcohol LET OP! De reflector- folie vereist andere reinigingsmiddelen dan het segment zelf.	Beschadiging. Zie paragraaf <b>"Fluent ID-reflectorfolie ver-</b> vangen" [▶ 138].	
Runners	Reinigen	DI-water, alcohol, mild reinigingsmiddel, ontsmettingsmiddel, oppervlaktereinigings- middel, mild zuur, base, bleek, DNAzap	Zie paragraaf <b>"Runners and segmenten reinigen"</b> [▶ 136].	
Vaste naaldenblok (MCA)	Parkeren, reinigen, controleren en afdek- ken	Alcohol	Zie paragraaf "Vastenaal- denblok (MCA 384) reinigen" [▶ 151].	
			VOORZICHTIG! Tecan advi- seert om het naaldenblok na elke ploeg te parkeren en van het instrument te verwijderen. Zorg altijd voor een correcte reiniging en correcte opslag van het naaldenblok. Controleer op verplaatsing van de naald.	



Instrument/ component	Systeemverzorgings- taak	Reinigingsmiddel/ wegwerpproduct/ apparaat	Referentie/ systeemverzorgingsactivi- teiten
Vaste naalden	Reinigen	Alcohol, bleekmiddel, pluisvrije doek	Zie paragraaf "Vaste punten reinigen" [▶ 131].
Wegwerpnaaldkegels	Reinigen	Alcohol, pluisvrije doek	Zie paragraaf "Wegwerp- punt-kegel reinigen" [▶ 130].
Was- en afvoerstation (Liquid FCA)	Reinigen	Dl-water, alcohol, mild reinigingsmiddel, ontsmettingsmiddel	Zie paragraaf "Wegwerp- punt-afvoer en wasstation- eenheid reinigen" [> 140].
Wegwerpnaald-af- voerlade en afvoeraf- dekkingen	Reinigen	DI-water, alcohol, mild reinigingsmiddel, ontsmettingsmiddel	Zie paragraaf "Wegwerp- punt-afvoerlade reinigen" [▶ 141].
Vloeistofsysteem (Liquid FCA)	Spoelen	Systeemvloeistof LET OP! Als het vloei- stofsysteem een hoog zoutgehalte heeft, spoelen met gedeïo- niseerd water.	Voer de methode Liquid FCA Routine Flush Main- tenance (Routinematig doorspoelonderhoud FCA- vloeistof) uit.
MCA-wasstation (MCA met vaste naal- den)	Reinigen	Alcohol, bleekmiddel, Decon 90, Contrad LET OP! Oppervlak- tereinigingsmiddelen, zoals Decon of Con- trad kunnen het pro- ces beïnvloeden. Daarom moet het proces bij gebruik van dergelijke reinigings- middelen zorgvuldig worden gevalideerd.	Voer de opdracht <b>Prime</b> <b>Wash Station (MCA384)</b> (Wasstation primen) uit.



Instrument/ component	Systeemverzorgings- taak	Reinigingsmiddel/ wegwerpproduct/ apparaat	Referentie/ systeemverzorgingsactivi- teiten
Wegwerpnaald-af- voerzak	Vervangen	Aanbevolen zakspeci- ficaties: B x L: 300 mm x 600 mm; dikte: 0,5 mm Materiaal: polypropy- leen, polyethyleen of copolymeer (autocla- veerbaar) LET OP! De afvoerzak moet voldoen aan de plaatselijke veilig- heidsrichtlijnen.	Zie paragraaf "Wegwerp- punt-afvoerlade reinigen" [▶ 141].
Systeemvloeistofre- servoir (Liquid FCA)	Ervoor zorgen dat de- ze schoon is	Systeemvloeistof	Zie paragraaf "Het systeem- vloeistofreservoir en afvoerre- servoir aansluiten" [> 146].
Afvoerreservoir (Liquid FCA en MCA met vaste naalden)	Legen en reinigen	DI-water, alcohol, mild reinigingsmiddel, oppervlaktereinigings- middel, desinfectie- middel, base, bleek	Afhankelijk van uw lokale laboratoriumregelgeving/ wetgeving dagelijks of we- kelijks reinigen. Zie paragraaf <b>"Het systeem-</b> vloeistofreservoir en afvoerre- servoir aansluiten" [> 146].
Veiligheidspaneel	Reinigen	DI-water, alcohol, mild reinigingsmiddel	Zie paragraaf "Veiligheidspa- nelen reinigen" [▶ 140].
Buisrotator	Oppervlakken, neer- houder en wasstation reinigen	Pluisvrije doeken met 2% bleekmiddel, 70% ethanol of 100% isopropanol	Zie paragraaf "De buisrotator reinigen" [▶ 131].
Doorpriknaald	Controleer de door- priknaalden visueel op eventuele afzettingen. Indien nodig reinigen.	70% ethanol of 2% bleekmiddel en pluisvrije doek	Zie paragraaf "Doorprikpun- ten reinigen" [▶ 131].
Doorpriknaald was- station en afvalslang	Reinig het wasstation en de slangen van het resterende monster- materiaal.	DI-water, mild reini- gingsmiddel, ont- smettingsmiddel. Gebruik geen bleek- middel zonder de on- derdelen van het wasstation na te spoelen met water.	Voer een uitgebreide spoe- ling van het wasstation uit, inclusief alle compartimen- ten (voor, midden en ach- terkant van het wasstation).



Instrument/ component	Systeemverzorgings- taak	Reinigingsmiddel/ wegwerpproduct/ apparaat	Referentie/ systeemverzorgingsactivi- teiten
Frida Reader	Verwijder het inzetstuk en bevestig de blinde stop in de Frida Rea- der	_	Zie paragraaf <b>"Frida Reader"</b> [▶ 153].

### 7.4.2 Wekelijkse systeemverzorging

Wekelijkse systeemverzorging moet op de laatste werkdag van elke week worden uitgevoerd.

Voer de methode **WeeklySystemCare** uit, indien beschikbaar gesteld door de keyoperator, of voer, naast de dagelijkse taken, elke individuele taak uit die van toepassing is op uw Fluent-armconfiguratie, in chronologische volgorde zoals vermeld in de onderstaande tabel.

Tab	21.	Tobol	voor	wokali	ikaa (	ovotoom	vorzoraina
i av.	24.	Iaper	VUUI	WEVEI	jnse :	sysieem	verzorging

Instrument/ component	Systeemverzorgings- taak	Reinigingsmiddel/ wegwerpproduct/ apparaat	Referentie/ systeemverzorgingsactivi- teiten
Wassysteem (MCA 384 met vaste naalden)	Buis en filter visueel controleren op vuil	Filter	Vuil filter. Raadpleeg de pa- ragraaf "Klantsupport" [▶ 203].
Pipetteerkop (MCA 384)	teerkop Een lektest uitvoeren -		Voer de methode MCA 384 leakage and zero dispense uit.
Plaatadapter (MCA 384)	Reinigen	Alcohol Perslucht (voor drogen)	Zie paragraaf " <b>Plaatadapter</b> reinigen (MCA 384)" [▶ 151].
Vloeistofsysteem (Liquid FCA)	Reinigen	Afhankelijk van de door de Fluent ver- werkte vloeistof Decon, Contrad, ba- se, mild zuur, desin- fectiemiddel Gevolgd door spoelen met water, alcohol en systeemvloeistof	Zie paragraaf "Vloeistofpad reinigen" [▶ 145].



Instrument/	Systeemverzorgings-	Reinigingsmiddel/	Referentie/
component	taak	wegwerpproduct/ apparaat	systeemverzorgingsactivi- teiten
Liquid FCA	Controleer de correcte bevestiging van injec- tiespuiten aan de kle- pinterface, en de cor- recte bevestiging van de injectiespuitplunjer bij de vergrendelings- schroef van de plun- jer.	_	Zie paragraaf <b>"Bevestiging</b> van injectiespuiten controle- ren" [▶ 146]
DiTi-kegels	Bevestiging van DiTi- kegel controleren	_	Zie paragraaf <b>"Een DiTi cone</b> vastzetten" [▶ 152]
Liquid FCA	Een lektest uitvoeren (Liquid FCA)	-	Voer de methode <b>Liquid</b> FCA-lektestuit.
Air FCA	Een lektest uitvoeren (Air FCA)	_	Voer de methode <b>Air FCA-</b> lektest en cLLD-zelftestuit.
Air FCA MultiSense	De MultiSense Air FCA cLLD-zelftest uitvoeren	-	Voer de methode <b>Air FCA-</b> lektest en cLLD-zelftestuit.
Systeemvloeistofre- servoir	Reinigen	DI-water, alcohol, mild reinigingsmiddel, oppervlaktereinigings- middel, desinfectie- middel, base, bleek	Zie paragraaf "Het systeem- vloeistofreservoir en afvoerre- servoir reinigen" [> 146].
Wasstation (Liquid FCA)	Reinigen	Reinigingsmiddel of antiseptische oplos- sing	-
RGA-grijpervinger- pads	Verwijder deeltjes en resten van grijpervin- gerpads	Pluisvrije doek met alcohol	Afvegen met reinigingsmid- del.
Dokstation en grijper- vingers (bevestigings- interface)	Verwijder deeltjes en resten van grijpervin- ger-bevestigingsinter- face (PCBA, magneet en conus)	Pluisvrije doek met alcohol	Afvegen met reinigingsmid- del.
Standalone-barco- descannervenster	Reinigen	Mild reinigingsmiddel	WAARSCHUWING! Kijk niet in de laserstraal. Zie de handleiding van de fabrikant van de barco- descanner. Zie paragraaf "Laserstraling instrument" [> 38].



Instrument/ component	Systeemverzorgings- taak	Reinigingsmiddel/ wegwerpproduct/ apparaat	Referentie/ systeemverzorgingsactivi- teiten
Scannervenster Flu- ent-ID en buisrotator	Op vuil en beschadi- ging controleren Indien nodig reinigen	Mild reinigingsmiddel DI-water om te spoe- len	WAARSCHUWING! Kijk niet in de laserstraal. Reinigen en spoelen met een zachte doek.
Reflector Fluent-ID en buisrotator	Op vuil en beschadi- ging controleren Indien nodig reinigen Mild reinigingsmiddel DI-water om te spoe- len		WAARSCHUWING! Kijk niet in de laserstraal. Reinigen en spoelen met een zachte doek.
FCA-grijper	Reinigen	Alcohol	-
Buisrotator	Aanwezigheid en be- vestiging van vergren- del- en positioneer- pennen controleren. Indien nodig, de pen- nen bevestigen of vervangen	_	Zie paragraaf "Vergrendel- pennen en positioneerpennen vervangen" [▶ 198].
MCA 96	Een lektest uitvoeren	_	Voer de methode <b>MCA 96-</b> <b>lekkage-</b> methode uit
MCA 96	Voer een pipetteer- prestatietest uit	-	<ul> <li>Voer de methode MCA 96 pipetteerprestatie- methode uit als:</li> <li>de levensduur van de kegelafdichting 90% is</li> <li>er naalden van 10 µl of 50 µl worden gebruikt</li> <li>Fluent Control geeft een waarschuwing zodra 90% van de levensduur van de kegelafdichting is bereikt.</li> </ul>

### 7.4.3 Maandelijkse systeemverzorging

De volgende tabel somt in chronologische volgorde de maandelijkse systeemverzorgingstaken op:



#### Tab. 25: Tabel voor maandelijkse systeemverzorging

Instrument/ component	Systeemverzorgings- taak	Reinigingsmiddel/ wegwerpproduct/ apparaat	Referentie/ systeemverzorgingsactivi- teiten
Software	De computer herstar- ten	_	De computer uitschakelen. Wacht 10 seconden. De computer weer inschakelen.
Armgeleiding	Reinigen	Wattenstaafje of een pluisvrije doek op een schroevendraaier	Zie paragraaf "Armgeleiding reinigen" [> 152].
MCA 96	Reinig MCH 96	Pluisvrije doek met alcohol, perslucht	Zie paragraaf " <b>Reinig MCH</b> 96" [▶ 154]

### 7.4.4 Regelmatige systeemverzorging



De intervallen waarop deze taken moeten worden uitgevoerd, moeten door de keyoperator worden bepaald.

De volgende tabel somt in chronologische volgorde de systeemverzorgingstaken op:

Tab. 26: Tabel voor periodieke systeemverzorging

Instrument/ component	Systeemverzorgings- taak	Reinigingsmiddel/ wegwerpproduct/ apparaat	Referentie/ systeemverzorgingsactivi- teiten
Kegel-hulsverbinding	Deeltjes verwijderen Oppervlakken reinigen	Alcohol, pluisvrije doek	
UVC-licht	Controleren op vin- gerafdrukken. Indien nodig reinigen.	Alcohol, pluisvrije doek	
MCA 96-grijpervingers	Visuele inspectie van grijpervingers, met name na botsingen van grijpervingers	Controleer de grijper- vingers op vuil en be- schadigingen. Als ze vuil zijn, maakt u ze schoon met alcohol en een pluisvrije doek.	

### 7.4.5 Jaarlijkse systeemverzorging

De jaarlijkse systeemverzorging helpt de nauwkeurigheid en precisie te behouden en de stilstandtijden van het instrument te minimaliseren. De verzorging ondersteunt ook een langere levensduur van de Fluent.



Neem contact op met de plaatselijke Tecan-serviceorganisatie om een vaste, jaarlijkse systeemverzorging af te spreken. Raadpleeg de paragraaf "Klantsupport" [> 203].

### 7.4.6 Tweejaarlijkse systeemverzorging

De volgende preventieve onderhoudstaken moeten elke 2 jaar worden uitgevoerd:

Tab.	27:	Tweeja	arlijkse	systeem	/erzorging
		,	,	,	55

component	Taak	Verwijzing
FCA-grijper	FCA-grijpervingers vervan- gen. Reset de teller in Fluent- Control.	Voor de bestelgegevens, zie de referentiehandleiding. Zie "Referentiedocumenten" [▶ 9].

### 7.5 Systeemverzorgingsactiviteiten

Om de systeemverzorgingstaken zoals beschreven uit te voeren, als volgt te werk gaan:

- Wissel naar systeemverzorgingsmodus. Zie paragraaf "Systeemverzorgingsmodus" [> 115].
- Neem de instructies zoals hieronder beschreven in acht.

### 7.5.1 Het instrument op een kabinet binnen het laboratorium verplaatsen

### **A VOORZICHTIG**

### Beschadiging van het kabinet!

Kabinetlegborden kunnen zijn verwijderd, bijvoorbeeld voor een centrifuge-installatie of afvalbak. Door verplaatsen van het instrument dat op een kabinet is geplaatst zonder geïnstalleerde legborden kan het kabinet worden beschadigd en letsel veroorzaken.

- Installeer de kabinetlegborden, alvorens het instrument te verplaatsen.
- Verplaats de kast alleen op een vlakke vloer zonder treden of sporen. Als er treden of sporen worden aangetroffen, gebruik dan de Fluent-hefstangen om het systeem over de obstructie te tillen of neem contact op met uw servicevertegenwoordiger.

Verplaats het instrument als volgt op een kabinet binnen een ruimte:

1. Zorg ervoor dat het kabinet veilig is geparkeerd en tegen wegrollen is beveiligd.



2. Zorg ervoor dat de kabinetlegborden (A) zijn geïnstalleerd.



- 3. Draai de moer van de kabinetpoot (B) met een steeksleutel.
- 4. Draai de rode schroef (D) op de kabinetpoten (C) om de vergrendeling te ontgrendelen en de wielen kunnen bewegen.



- 5. Verplaats het instrument op het kabinet naar de nieuwe locatie.
- 6. Zorg ervoor dat het kabinet veilig is geparkeerd en tegen wegrollen is beveiligd.

### 7.5.1.1 Het instrument waterpas zetten

Zet het instrument als volgt waterpas:





1. Laat met een steeksleutel alle verstelbare poten neer tot de kabinetwielen handmatig kunnen worden gedraaid.

2. Maak de borgmoer (A) van de bijbehorende voet los.



3. Plaats het referentiesegment overeenkomstig de hieronder aangegeven rasterposities.

Instrumentmaat 480: Rasterpositie links 1 en rasterpositie rechts 21. Instrumentmaat 780: Rasterpositie links 1 en rasterpositie rechts 41.





Instrumentmaat 1080: Rasterpositie links 1 en rasterpositie rechts 59.

- 4. Gebruik de waterpas om ervoor te zorgen dat het instrument horizontaal en verticaal is uitgelijnd.
- 5. Stel het kabinetniveau af zoals vereist (rechtsom omhoog, linksom omlaag).



- 6. Na het waterpas maken van het instrument, de borgmoeren op de kabinetpoten weer vastdraaien.
- 7. Zorg ervoor dat het kabinet veilig is geparkeerd en tegen wegrollen is beveiligd.

### 7.5.2 Wegwerppunt-kegel reinigen

Reinig de wegwerppunt-kegel als volgt:

- 1. Reinig de wegwerppunt-kegels met alcohol met een pluisvrije doek.
- Controleer de wegwerppunt-kegels en uitstekende punt tijdens systeemverzorging.
   Voor Liquid ECA: Controleer of de buisverlenging die uit de buitenz

Voor Liquid FCA: Controleer of de buisverlenging die uit de buitenzijde van de kegel uitsteekt niet is beschadigd.

3. Zorg ervoor dat buisverlengingen schoon en vrij van afzettingen zijn.



### 7.5.3 Vaste punten reinigen

### **⚠ VOORZICHTIG**

### Risico op letsel door vaste punten tijdens de reiniging

Pipetteren met vaste punten kan letsel veroorzaken.

• Contact met pipetteerpunten en aerosolen bij toegang tot de werktafel vermijden door geschikte beschermende kleding te dragen.

Reinig de vaste punten als volgt:

- 1. Reinig de vaste punten met alcohol en een pluisvrije doek.
- 2. Zorg ervoor dat vaste punten schoon en vrij van afzettingen zijn.

### 7.5.4 Doorprikpunten reinigen

Reinig de doorprikpunten met de **doorprikpunt-reinigingsonderhoud** methode. Deze methode moet aan uw werktafelindeling worden aangepast.

Het script omvat de volgende stappen:

- 1. De werktafel voorbereiden (bijv. laboratoriummateriaal en hardware).
- 2. Doorprikken tot Z-start van 8 lege buizen met dop op een buisrotator of op een buisneerhouder-drager.
- 3. Reinig het toegankelijke deel van de doorprikpunten handmatig met 70% ethanol of 2% bleek op een pluisvrije doek. Voorkom contact met de scherpe punt van de doorprikpunten.
- 4. Voer wastaken na handmatige reiniging uit.

### 7.5.5 De buisrotator reinigen

### Algemene reinigingsprocedure

- 1. Reinig alle delen van de buisrotator met pluisvrije doeken en dompel ze in een van de volgende reinigingsvloeistoffen: 2% bleekmiddel, 70% ethanol, 100% isopropanol
- 2. Veeg de onderdelen schoon met de gedrenkte doeken om ze te reinigen en te ontsmetten.

Gebruik wattenstaafjes voor delen die niet bereikbaar zijn met een pluisvrije doek.

3. Veeg de reinigingsvloeistoffen binnen 5 minuten na het aanbrengen weg met in water gedrenkte doeken.



### Buis-neerhouderplaat verwijderen en reinigen

1. Om de neerhouderplaat (A) los te maken, de plaat met een hand vasthouden en de neerhouder-vergrendelpen (B) met de andere hand omlaag trekken.



2. Verwijder de neerhouderplaat van de buisrotator.



3. Reinig de neerhouderplaat volgens de algemene, bovenstaande instructies. Als alternatief kan de buisneerhouder worden geïncubeerd in een bad met 2% bleek gedurende maximaal 2 uur.

#### De buisrotatoroppervlakken reinigen

1. Reinig de toegankelijke oppervlakken van de buisrotator volgens de algemene, bovenstaande instructies.





2. Om de trommelpositie handmatig te wijzigen, de trommel met een hand vasthouden en de magneetspoel-vrijgavetoets indrukken.

- 3. Roteer de trommel handmatig en laat de magneetspoel-vrijgavetoets weer los.
- 4. Roteer de trommel tot deze door de magneetspoel wordt vergrendeld.
- 5. Reinig de oppervlakken die eerder volgens de algemene, bovenstaande instructies niet toegankelijk waren.

### De neerhouderplaat monteren

- 1. Plaats de neerhouderplaat boven op de buisrotatortrommel.
- 2. Druk met de ene hand de neerhouderplaat naar de onderzijde van het instrument en duw de zwarte schuif naar achteren om de neerhouderplaat in positie te vergrendelen.





### 7.5.6 Het buisrotator-wasstation reinigen

#### Algemene reinigingsprocedure

- Het wasstation kan worden gereinigd op de werktafel of kan voor reiniging worden gedemonteerd.
- ✓ Gebruik voor betere toegang in plaats van een doek een flessenborstel.
- 1. Reinig alle delen van het wasstation met pluisvrije doeken en dompel ze in een van de volgende reinigingsvloeistoffen: 2% bleekmiddel, 70% ethanol, 100% isopropanol
- 2. Reinig en ontsmet de onderdelen met gedrenkte doeken.
- 3. Veeg de reinigingsvloeistoffen binnen 5 minuten na het aanbrengen weg met in water gedrenkte doeken.

### Buisrotator-wasstation demonteren

1. Duw het wasstation-vrijgavehendel richting de barcodescannerbehuizing, en til het wasstation met de andere hand op.







2. Ontkoppel de afvoerbuis en plaats de connectors in de afvoerbuishouders.

#### **Buisrotator-wasstation monteren**

1. Sluit de afvoerbuisconnectors aan.







 Monteer het wasstation op de geleide-assen (A) en druk het op de basisplaat. Controleer of de vrijzethendel op zijn plaats terugspringt en het wasstation in positie houdt.

### 7.5.7 Runners and segmenten reinigen

Reinig runners en segmenten als volgt:

### LET OP

### Storing van vloeiststofdetectie (cLLD)!

Mogelijke storing van de vloeiststofdetectie (cLLD) door verminderd contact tussen loper en deksegment.

Zorg er altijd voor dat de runners en segmenten schoon en droog zijn.

1. Verwijder de runners van het instrumentdek.

Segmenten en nesten ter plaatse gereinigd.

2. Wis het oppervlak van de runners, segmenten en nesten schoon met reinigingsmiddel.

Spoel de runners, segmenten en nesten schoon met DI-water.

3. Plaats de runners terug op het instrumentdek.

### 7.5.8 Dektrays reinigen

Reinig de dektrays als volgt:

- ✓ Segmenten boven dektray worden verwijderd. Zie "Segmenten verwijderen" [▶ 85].
- ✓ Als deksegmenten, zoals Fluent ID of MCA Active Carrier niet kunnen worden verwijderd, de dektrays op een open dekpositie schuiven.
- 1. Verwijder de dektrays van het instrument.



- 2. Leeg de trays door verwijderen van de vloeistof volgens het laboratoriumprotocol voor deze specifieke vloeistof.
- 3. Als de dektrays worden beschadigd of verloren gaan, moeten ze worden vervangen.
- 4. Veeg het oppervlak van de dektrays schoon met het reinigingsmiddel.
- Plaats de dektrays terug in het instrument.
   Positioneer de dektrays zoals hieronder afgebeeld.
   Aangrenzende dektrays moeten in elkaar grijpen.



Afb. 44: Onjuiste dektraypositionering



Afb. 45: In elkaar grijpende dektrays





Afb. 46: Correcte dektraypositionering

### 7.5.9 Fluent ID-reflectorfolie vervangen

- ✓ Zelfhechtende reflectorfolie
- 1. Verwarm de reflectorfolie. Gebruik een verwarmingspistool.
- 2. Verwijder de reflectorfolie.



3. Verwijder alle resten met alcohol.





4. Breng nieuwe, zelfhechtende reflectorfolie aan op het bovenste einde van de reflector.

### 7.5.10 Fluent ID-reflectorfolie op DiTi-afvoerlade aanbrengen

- ✓ Zelfhechtende reflectorfolie
- 1. Breng tape aan op de DiTi-afvoerlade volgens onderstaande afbeelding.



2. Breng de zelfhechtende reflectorfolie aan op de DiTi-afvoerlade volgens onderstaande afbeelding.





De laserstraal moet zich in het centrum van de reflectorfolie bevinden.

3. Verwijder de tape van de DiTi-afvoerlade.

### 7.5.11 Veiligheidspanelen reinigen

Reinig de veiligheidspanelen als volgt:

1. Veeg het binnen- en buitenvlak van de veiligheidspanelen schoon met reinigingsmiddel.

### 7.5.12 Wegwerppunt-afvoer en wasstationeenheid reinigen

Reinig de wegwerppunt-afvoer en het wasstation als volgt:

- 1. Druk op de snelspantoets (B).
- 2. Schuif het wasstation naar achteren.



Afb. 47: Bevestiging voor het verwijderen van de zakbehuizing

3. Verwijder het wasstation van de wegwerppunt-afvoer en de wasstationeenheid.





Afb. 48: Het wasstation verwijderen

- 4. Veeg het oppervlak van het wasstation schoon met een reinigingsmiddel en verwijder gemorste reagentia volledig.
- 5. Druk op de snelspantoets (B).
- 6. Positioneer het wasstation (A).
- 7. Duw het wasstation vooruit.



Afb. 49: Breng het wasstation weer aan

### 7.5.13 Wegwerppunt-afvoerlade reinigen

Reinig de wegwerppunt-afvoerlade als volgt:

- ✓ Veiligheidspaneel aan de voorzijde is open.
- 1. Verwijder de afdekking (A) van de wegwerppunt-afvoerlade.





Afb. 50: De afdekking van wegwerppunt-afvoerlade verwijderen

2. Verwijder de wegwerppunt-afvoerlade (B) van de houder.



Afb. 51: Wegwerppunt-afvoerlade verwijderen

3. Houd een doek onder de opening aan de onderzijde van de wegwerppuntafvoerlade (C).

Voorkom druppelen van verontreinigde stoffen.



Afb. 52: Wegwerppunt-afvoerlade behandelen

- 4. Plaats de wegwerppunt-afvoerlade en de afdekking een met reinigingsmiddel gevulde bak.
- 5. Laat de lade 30 minuten tot 4 uur in de vloeistof.
- 6. Neem de wegwerppunt-afvoerlade en afdekking uit de bak en plaats ze op een droge doek.



- 7. Laat ze drogen.
- 8. Monteer de wegwerppunt-afvoerlade (B) weer op de houder.



- Afb. 53: Inzetstuk van wegwerppunt-afvoerlade weer monteren
- 9. Zorg ervoor dat de positioneerpen correct in de sleuf (D) is gestoken.
- 10. Plaats de afdekking (A) op de afvoerlade.



Afb. 54: Pen en afdekking positioneren

### 7.5.14 Wegwerppunt-afvoerzak vervangen

Reinig de wegwerpnaald-afvoerzak als volgt:

1. Til de bevestiging (A) op, en schuif zakbehuizing vooruit.



Afb. 55: Bevestiging voor het verwijderen van de zakbehuizing



- 2. Verwijder de wegwerpzakbehuizing (A).
- 3. Verwijder de wegwerpnaald-afvoerzak (B).
- 4. Voer de wegwerpnaald-afvoerzak af volgens de laboratoriumrichtlijnen.
- 5. Plaats de nieuwe wegwerpzak (B) in de lege zakbehuizing (B).



Afb. 56: Zakbehuizing en wegwerpnaald-afvoerzak



Afb. 57: Juiste montage van de afvalkoker op de werktafel



### 

Een verkeerd geplaatste afvalkoker kan ervoor zorgen dat de arm botst en/of de naalden verkeerd worden uitgeworpen. Zorg ervoor dat de afvalkokers correct zijn geplaatst zoals hieronder weergegeven: voeg afbeeldingen toe die aangeven of de afvalkokers correct/onjuist zijn geplaatst

6. Schuif de zakbehuizing in de juiste positie, sluit hem dan met bevestiging (A).




Afb. 58: Bevestiging sluiten

#### 7.5.15 Vloeistofpad reinigen

- Verwijder proteïneresten van de binnenzijde van vaste naalden met mild zuur en daarna een basisch reinigingsmiddel.
- Verwijder nucleïnezuurresten van de binnenzijde van vaste naalden met een basisch reinigingsmiddel.
- Reinigingsmiddelen, zoals Decon/Contrad, kunnen het proces beïnvloeden. Daarom moet het proces bij gebruik van dergelijke reinigingsmiddelen zorgvuldig worden gevalideerd.
- Isopropanol is een uiterst effectief ontsmettingsmiddel. Het verdampt snel. Daarna zijn de oppervlakken onmiddellijk klaar voor gebruik.
- Gebruik alleen goedgekeurde reinigingsmiddelen. Gebruik geen bleekmiddeloplossingen om het gehele vloeistofsysteem te spoelen.

Reinig het vloeistofpad als volgt:

- 1. Ontkoppel de systeemvloeistofbuis van het vloeistofreservoir.
- 2. Sluit de onderhoudsbuis (30043739) aan op de systeembuis.
- 3. Plaats het open einde van de onderhoudsbuis in een fles met reinigingsmiddel.
- 4. Spoel met reinigingsmiddel (20 ml met RapidWash en 10 ml met verdunner).
- 5. 20 minuten in de vloeistof laten zitten.
- 6. Plaats de buis in een fles met DI-water.
- 7. Spoel twee keer met DI-water (20 ml met RapidWash en 10 ml met verdunner).

## 

#### Ontvlambare vloeistoffen!

Brandgevaar door ontvlambare vloeistoffen of systeemvloeistof.

- Vermijd de vorming en ophoping van brandbare dampen.
- Bedien het systeem niet zonder dektrays.
- 8. Plaats de buis in een fles met alcohol.
- 9. Spoel met alcohol (20 ml met RapidWash en 10 ml met verdunner).
- 10. Verwijder de onderhoudsbuis uit de systeembuis en sluit de systeembuis aan op het systeemvloeistofreservoir.



- 11. Spoel twee keer met DI-water (20 ml met RapidWash en 5 keer met het verdunnervolume).
- 12. Controleer op bellen in de buis.
- 13. Spoel opnieuw, als er bellen zichtbaar zijn.

#### 7.5.16 Het systeemvloeistofreservoir en afvoerreservoir aansluiten

Bereid het systeemvloeistofreservoir en het afvoerreservoir als volgt voor:



Storingsvrij bedrijf is alleen gegarandeerd, als de originele reservoirs met het besturingssysteem van Tecan worden gebruikt. Voor het eerste gebruik moet het systeemvloeistofreservoir handmatig grondig worden schoongespoeld om eventuele vaste deeltjes uit de fles te verwijderen. Zie paragraaf "Het systeemvloeistofreservoir en afvoerreservoir reinigen" [> 146].

- ✓ Tecan-reservoir met een capaciteit van meer dan 20 liter
- 1. Zorg ervoor dat het vloeiststofdetectiesysteem (A, B) correct is aangesloten.
- 2. Zorg ervoor dat de buizen (C, D) correct zijn aangesloten.



#### 7.5.17 Het systeemvloeistofreservoir en afvoerreservoir reinigen

Reinig het vloeistofreservoir en het afvoerreservoir als volgt:

- 1. Leeg het wasvloeistofreservoir handmatig.
- 2. Reinig het vloeistofreservoir in een bak met reinigingsmiddel, en spoel het dan schoon.
- 3. Ontsmet het vloeistofreservoir met alcohol.
- 4. Sluit het systeemvloeistof- en het afvoerreservoir aan, zie paragraaf "Het systeemvloeistofreservoir en afvoerreservoir aansluiten" [▶ 146].

#### 7.5.18 Bevestiging van injectiespuiten controleren

Controleer de correcte bevestiging van injectiespuiten als volgt:

#### 7 - Systeemverzorging Systeemverzorgingsactiviteiten





Afb. 59: Bevestiging controleren

A Zuiger-vergrendelingsschroef B Injectiespuitschroef

- Verplaats de plunjers naar het midden van de spuiten met behulp van een onderhoudsscript dat lucht opzuigt.
   **Opmerking:** Het onderhoudsscript moet worden geleverd door de FluentControl-beheerder van het laboratorium.
- 2. Draai de injectiespuitschroef (B) aan dat wil zeggen rechtsom draaien.
- 3. Draai de zuiger-vergrendelschroef (A) rechtsom om hem vast te draaien.

## 7.5.19 Pakkingen controleren (MCA 384)

Controleer pakkingen (MCA) als volgt:

1. Controleer de pakkingen (A) op beschadiging.



Controleer of de pakkingen niet beschadigd zijn. Een beschadigde pakking moet worden vervangen. Zie paragraaf "Pakkingen vervangen (MCA 384)" [▶ 148].



## 7.5.20 Pakkingen vervangen (MCA 384)



Afb. 60: Onderdelen en gereedschappen

- A Pakking
- C Stompe buis-insteekgereedschap
- E Stompe buis-verwijderingsgereedschap
- B Stompe buis

D

- Pakking-insteekgereedschap
- Vervang pakkingen (MCA) als volgt:
- 1. Verwijder de adapterplaat.
- 2. Beweeg de pipetteerkop naar voren en til hem zo hoog mogelijk op.
- 3. Duw het stompe buis-verwijderingsgereedschap ca. 2 mm in het kanaal.





4. Borg de stompe buis door het gereedschap rechtsom te draaien.

5. Trek de stompe buis uit het kanaal.







6. Verwijder de pakking met een langbektang uit het kanaal.

- 7. Smeer de nieuwe pakking met een film van minerale olie.
- 8. Steek de gesmeerde pakking met het pakking-insteekgereedschap in het kanaal.



9. Behandel het buitenoppervlak van een nieuwe stompe buis met minerale olie.



10. Steek de gesmeerde stompe buis met het stompe buis-insteekgereedschap in het kanaal.

De stompe buis moet volledig in het kanaal gestoken zijn.



- 11. Controleer de nieuwe pakking om uitlijning met de andere pakkingen te garanderen.
- 12. Voer de **MCA-lektest** methode.

#### 7.5.21 Vastenaaldenblok (MCA 384) reinigen

- 1. Reinig het vastenaaldenblok (MCA) met alcohol en een pluisvrije doek.
- 2. Veeg het naaldenblok droog met een pluisvrije doek of blaas het naaldenblok droog met olievrije perslucht.
- 3. Zorg ervoor dat het vastenaaldenblok (MCA) schoon en vrij van afzettingen is.
- 4. Bewaar het naaldenblok in de naaldenblokdoos.

Om besmetting bij het werken met het naaldenblok te voorkomen:

- Het naaldenblok moet op een stofvrije plaats worden bewaard.
- Raak de naalden nooit aan met uw vingers. Het naaldenblok moet tijdens het werk altijd door het PEEK-blok worden vastgehouden.
- Plaats het naaldenblok nooit met de naalden op de tafel.

## 7.5.22 Plaatadapter reinigen (MCA 384)

Reinig de plaatadapter als volgt:

- 1. Controleer op vuil in de kanalen.
- 2. Verwijder vet met alcohol.
- 3. Droog met perslucht.



## 7.5.23 Armgeleiding reinigen

Reinig de armgeleiding als volgt:

- 1. Reinig de armgeleidingsroller (A) op de armgeleiding met een wattenstaafje of een pluisvrije doek op een schroevendraaier.
- 2. Reinig de armrails (B) met een pluisvrije doek.
- 3. Reinig het bovenste oppervlak van de geleiderail op de MCA-armgeleiding met een pluisvrije doek.



## 7.5.24 Een DiTi cone vastzetten

Zet de Air FCA DiTi cone als volgt vast:

- 1. Houd de puntadapter (D) en de punt-uitwerpbuis (C) vast.
- 2. Draai de DiTi cone (A) aan met de DiTi-steeksleutel (B).



3. Voer de FCA routineonderhoud methode.



### 7.5.25 Frida Reader

Inzetstuk



*Afb. 61:* Frida Reader-inzetstuk Plaats het inzetstuk voor montage in de Frida Reader en lijn de markeringen uit.

#### Blinde stop



Afb. 62: Frida Reader-blinde stop

De blinde stop beschermt de Frida Reader, als het inzetstuk is verwijderd. Bevestig de blinde stop voor montage in de Frida Reader.



## 7.5.26 Reinig MCH 96

Controleer de uitwerpplaat (A) op zichtbare schade. Morsen van bepaalde procesvloeistoffen (zoals DMSO of acetonitril) of reinigingsmiddelen (zoals bleekmiddel) kan leiden tot beschadiging van de plaat. Als dit het geval is, neem dan contact op met uw lokale serviceorganisatie voor vervanging.



- 1. Gebruik het verplaatsingsgereedschap om de uitwerpas naar beneden te drijven. 2. . 3.
- 2. Gebruik olievrije perslucht om de bovenkant van de uitwerpplaat (C), de kegelplaat en de kegels (B) te reinigen
- 3. Gebruik een pluisvrije doek en alcohol om de onderkant van de uitwerpplaat (D) te reinigen.





De linkerafbeelding hieronder toont de uitwerpplaat vóór het reinigen en de rechterafbeelding toont de uitwerpplaat na het reinigen.







# 8 Verhelpen van storingen

Lees dit hoofdstuk voor informatie over het voortzetten van de bediening na een probleem met de Fluent. Voor meer informatie of in het geval van problemen die niet in deze handleiding worden behandeld of die onvoldoende gedetailleerd zijn, zie de paragraaf "Klantsupport" [> 203].

# 8.1 Veiligheidsinstructies voor dit hoofdstuk

# **▲ VOORZICHTIG**

## Kruisbesmetting door beschadigde punten na botsen!

Gebogen punten of beschadigde puntcoating veroorzaken onnauwkeurigheid bij het pipetteren en fouten bij vloeistofdetectie.

 Controleer de vaste punten na een botsing. Zie paragraaf "Vaste punten controleren" [▶ 179].

# 8.2 Tabellen voor het verhelpen van storingen

De tabellen voor het verhelpen van storingen sommen mogelijke problemen, oorzaken en correctiemaatregelen op. Voor meer informatie of in het geval van problemen die niet in deze handleiding worden behandeld of die onvoldoende gedetailleerd zijn, zie de paragraaf "Klantsupport" [> 203].

## 8.2.1 Verhelpen van storingen van het instrument

Tab. 28: Tabel voor het verhelpen van storingen van het instrument

Probleem/fout	Mogelijke oorzaak	Correctiemaatregel
Lekkage van systeemvloei- stof	Buis en/of buisverbin- dingen zijn niet cor- rect bevestigd. Injectiespuit lekt.	Raadpleeg de paragraaf <b>"Klantsupport</b> " [▶ 203].
Communicatie- fout	Voeding is niet AAN. Voeding of communi- catie is onderbroken. Geen communicatie.	Schakel het instrument uit. Wacht tot de instrument-statuslamp en stroomvoorzieningslamp uitgaan. Schakel de pc uit. Controleer kabel en stekkers. Schakel het instrument en de pc in.
	X-, Y- of Z-aandrij- ving geblokkeerd.	Controleer op hindernissen. LET OP! Zorg ervoor dat de armen vrij kunnen bewegen.



Probleem/fout	Mogelijke oorzaak	Correctiemaatregel
Initialisatiefout	Armen kunnen niet worden geïnitialiseerd.	Controleer op hindernissen. LET OP! Zorg ervoor dat de armen vrij kunnen bewegen.
	Hardwareprobleem.	Raadpleeg de paragraaf "Klantsupport" [> 203].
Deursensor van veiligheidspa- neel aan de voorzijde en deurvergrende- ling zijn be- schadigd	Mechanische fout van de deurvergrendelin- gen.	Schakel het instrument uit. Raadpleeg de paragraaf <b>"Klantsupport"</b> [▶ 203].
Veiligheidspa- neel ontbreekt of is bescha- digd	Veiligheid kan niet worden gegaran- deerd.	Schakel het instrument uit. Raadpleeg de paragraaf "Klantsupport" [▶ 203].
Vloeiststofde- tectie (cLLD) fout	Vuil contactoppervlak. Onvoldoende contact tussen laboratorium- materiaal en seg- ment.	Bereid het dek voor. Zie paragraaf . Reinig het contactoppervlak. Zie para- graaf "Runners and segmenten reinigen" [▶ 136].
	Systeemvloeistof heeft een geleidbaar- heid > 10 µS/cm voor cLLD-compati- biliteit.	Neem contact op met de key-operator.

## 8.2.2 Flexibele kanaalarm (FCA) verhelpen van storingen

Tab. 29: Tabel voor het verhelpen van storingen van de flexibele kanaalarm

Probleem/fout	Mogelijke oorzaak	Correctiemaatregel
Losse DiTi-ke- gel VOORZICH- TIG! Onnauw- keurige pipet- teervolumes!	Niet-correct bevestig- de DiTi-kegel.	Draai de DiTi-kegel aan.
Wegwerpnaald niet opgehaald	Niet-correct bevestig- de DiTi-kegel.	Draai de DiTi-kegel aan.



Probleem/fout	Mogelijke oorzaak	Correctiemaatregel	
Wegwerpnaald niet afgevoerd	Niet-correct bevestig- de DiTi-kegel.	Draai de DiTi-kegel aan.	
	Opnieuw gebruikte Di- Ti's	Zorg dat naalden nieuw zijn. Hergebruik van DiTi's wordt niet aanbevolen.	
		Wegwerpnaalden niet uitgeworpen in af- valkoker/afvalkoker is niet correct ge- plaatst	
Wegwerpnaal- den niet uitge- worpen in af- valkoker	De afvalkoker bevindt zich niet op de juiste plaats	Zorg ervoor dat de afvalkokers correct zijn geplaatst. Zie paragraaf "Wegwerp- punt-afvoerlade reinigen" [> 141]	
Naalden niet uitgelijnd ten opzichte van la-	Drager in verkeerde positie.	Zorg voor de correcte positie van de drager. Zie paragraaf <b>"Standaard runners</b> inladen" []> 87].	
boratoriumma- teriaal op een enkelvoudige drager	plaats vergrendeld. Laboratoriummateri- aal niet correct gepo- sitioneerd.	Vergrendel het segment op zijn plaats. Zie paragraaf " <b>Segment controleren</b> " [▶ 196].	
Naalden niet uitgelijnd ten opzichte van la- boratoriumma- teriaal op meerdere dra- gers	Defecte armuitlijning door een botsing.	Raadpleeg de paragraaf <b>"Klantsupport"</b> [ <b>▶ 203]</b> .	
Naald botst te- gen onderzijde van laboratori- ummateriaal	Onjuist laboratorium- materiaal. Laboratoriummateri- aal niet correct gepo- sitioneerd.	Zorg ervoor dat het laboratoriummateri- aal op het dek overeenkomt met de dekindeling van de methode.	
DiTi druppelt	Vuile DiTi-kegel ver- oorzaakt lekkage.	Reinig de DiTi-kegel.	
	Opnieuw gebruikte Di- Ti's	Zorg dat naalden nieuw zijn. Hergebruik van DiTi's wordt niet aanbevolen.	
Foutmelding: Druk buiten be- reik (Air FCA)	Nat inlinefilter na op- zuiging met onjuiste DiTi-maat.	Zorg ervoor dat de maat van de DiTi op het dek overeenkomt met de maat die in de methode is gedefinieerd. Zorg er- voor dat de Diti-kegels correct zijn vast- gedraaid Controleer het inlinefilter. Zie paragraaf	



Probleem/fout	Mogelijke oorzaak	Correctiemaatregel
Foutmeldingen:	Magnetisch veld	-
DiTi niet opge- haald	stoort DiTi-aanwezig- heidssensor.	
DiTi niet geval- Ien		
Vloeistofdetec- tie:	DiTi-kegel zit los	Zet de DiTi-kegel vast (zie <b>"Een DiTi co-</b> ne vastzetten" [▶ 152]).
naald detecteert geen vloeistof		
O-ringen ver- sleten	MultiSense-naal- dadapters	Vervang de O-ringen en X-ring op de MultiSense-naaldadapters.

## 8.2.3 Mengen en doorprikken

Tab. 30: Verhelpen van storingen

Symptoom	Mogelijke oorzaak	Correctiemaatregelen
Piercing tip kan met de software niet wor- den teruggetrokken	Piercing tip klemt	Zie paragraaf "Vastgeklemde doorprikpunten terugtrekken" [▶ 190].
Beschadigde piercing tip	Verbogen piercing tip Beschadigde punt	Vervang de piercing tip. Zie paragrafen "Doorprikpunten verwijderen" [▶ 183] en "Doorprikpunten installeren" [▶ 186].
Doorprikfouten	Piercing tip te droog	Smeren met water (wassta- tion)
	Beschadigde punt	Vervang de piercing tip. Zie
	Verbogen piercing tip	verwijderen" [▶ 183] en "Doorprikpunten installeren" [▶ 186].
	Onjuiste doorprikparameter	Neem contact op met de key-operator.
	Onjuist bewegingstype ge- bruikt	Neem contact op met de key-operator.
	Onjuiste buistypen gebruikt	Gebruik ondersteunde bui- zen. Zie paragraaf <b>"Buisrota-</b> tor-runners" [▶ 64].
	Arm heeft zijn levensduur bereikt	Raadpleeg de paragraaf "Klantsupport" [> 203].





	Symptoom	Mogelijke oorzaak	Correctiemaatregelen
	Problemen met vloei- stofverwerking	Verstopte piercing tips	Spoel piercing tips door. Controleer de wasprocedure in het algemeen.
		Beschadigde punt	Vervang de piercing tip. Zie paragrafen "Doorprikpunten verwijderen" [▶ 183] en "Doorprikpunten installeren" [▶ 186].
	Injectiespuiten niet correct gemonteerd	Controleer de correcte be- vestiging van de injectie- spuiten. Zie paragraaf <b>"Be-</b> vestiging van injectiespuiten controleren" [▶ 146].	
		Bellen in vloeistofsysteem	Spoelen. Raadpleeg de paragraaf "Klantsupport" [▶ 203].
	Hemolyseproblemen	Monsterverdunning	Hoger overtollig volume of afscheidingsvolume
			Zoutoplossing 0,9% als af- scheidingsvolume
			Lagere pipetteersnelheden
	Beschadigde punt	Vervang de piercing tip. Zie paragrafen "Doorprikpunten verwijderen" [> 183] en "Doorprikpunten installeren" [> 186].	
		Mengparameters	Ervoor zorgen dat de ge- bruikte rotatie/oscillatiepa- rameters van de buisrotator niet tot hemolyse leidt



Symptoom	Mogelijke oorzaak	Correctiemaatregelen
Monster in injectie- spuit	Willekeurig	Reinig het systeem. Zie pa- ragraaf <b>"Vloeistofpad reini-</b> gen" [▶ 145].
	Verkeerd luchtkussen.	Valideer wasprocedure.
	Injectiespuiten niet correct gemonteerd.	Controleer de correcte be- vestiging van de injectie- spuiten. Zie paragraaf <b>"Be-</b> vestiging van injectiespuiten controleren" [▶ 146].
		Voer de FCA-lekmethode.
		Groter eerste luchtkussen.
		Langzamere opzuigsnelhe- den.
	Onjuiste monstervoorberei- ding voor vloeistofverwer- king. Monsterbronbuizen bevatten vaste deeltjes zo-	Zorg voor een goede mon- stervoorbereiding om pipet- teren van monstervloeistof mogelijk te maken.
	ais stoiseis, ceiresten etc.	Zorg ervoor dat de mon- sterbronbuizen geen vaste deeltjes zoals stolsels, cel- resten etc. bevatten.
	Onjuiste monstervoorberei- ding voor vloeistofverwer- king. Buizen zijn niet goed gevuld en bevatten nog een	Zorg ervoor dat de mon- sterbronbuizen correct zijn gevuld met het doelvolume van de buis.
	eerste luchtkussen tijdens het doorprikken reduceert.	Controleer of de monster- bronbuizen geen vacuüm bevatten.
		Vergroot het eerste lucht- kussen om een eventueel resterend vacuüm te com- penseren.
Piercing tip verbuigt tijdens wasprocedure	Piercing tip is niet gecen- treerd in de reinigingsope- ningen van het wasstation	Maak een kopie van het wasstation en programmeer de pipetteerposities.



Symptoom	Mogelijke oorzaak	Correctiemaatregelen
Onjuist vloeistofni- veaudetectie: alleen op specifieke kanalen	Verbogen piercing tip: Pier- cing tip is verbogen raakt daarom de buiswand tijdens doorprikken	Vervang de piercing tip. Zie paragrafen "Doorprikpunten verwijderen" [▶ 183] en "Doorprikpunten installeren" [▶ 186].
	Doorprikpositie is onjuist, daarom raakt de piercing tip de buiswand tijdens door- prikken.	Gebruik laboratoriummateri- aal van Tecan. Zie para- graaf <b>"Buisrotator-runners"</b> [▶ 64].
		Pipetteerpositie van labora- toriummateriaal program- meren/afstellen
	Uitlijning van piercing tip is onjuist.	Monteer piercing tip met de opening richting de voorzij- de van het instrument. Zie paragraaf "Doorprikpunten installeren" [> 186].
Verkeerde vloeistofni- veaudetectie: con- stante afwijking van verwacht vloeistofni- veau en gedetecteerd vloeistofniveau	Productietoleranties van de arm, de buisrotator en/of het instrument in combinatie met de doorprikkracht kun- nen leiden tot een merkbare Z-offset tijdens vloeistofni- veaudetectie.	Teach/adjust the custom attribute "PiercingDetecti- onHeightCompensation" in the tube labware definition

## 8.2.4 Multikanaalarm (MCA) verhelpen van storingen

Tab. 31: Tabel voor het verhelpen van storingen van multikanaalarm

Probleem/fout	Mogelijke oorzaak	Correctiemaatregel
Punten zijn niet uitge- lijnd ten opzichte van de dragers	Mechanische fout	Raadpleeg de paragraaf <b>"Klantsupport"</b> [ <b>▶ 203]</b> .
	Armbotsing	
Multititerplaat en pi- petteerkop niet 100% parallel	Botsing	Raadpleeg de key-operator om de pa- rallelliteit van pipetteerkop en dek te controleren.
Tijdens het pipetteer- proces stopt de pi- petteerkop en gene- reert een fout Opzuig- en afgifteversnelling is snel vergeleken met de snelheid Opzuig- en afgifteverlangzamin is te snel vergeleken met snelheid	Opzuig- en afgifteversnelling is te snel vergeleken met de snelheid. Opzuig- en afgifteverlangzaming	Versnelling moet in een redelijke ver- houding staan tot de opzuig- en afgif- tesnelheid.
	is te snel vergeleken met snelheid.	Verlangzaming moet in een redelijke verhouding staan tot de opzuig- en af- giftesnelheid.
		Probleem kan niet worden verholpen. Raadpleeg de paragraaf <b>"Klantsupport"</b> [▶ 203].



Probleem/fout	Mogelijke oorzaak	Correctiemaatregel
Meerdere of alle pi- petteerkanalen lekken	Verkeerde wegwerppunten, vaste- puntblok, vastepuntadapter, punt- kegelafdichtingen of pakkingen.	Gebruik uitsluitend wegwerppunten, vastepuntblok, vastepuntadapter, punt- kegelafdichtingen of pakkingen van Te- can. Raadpleeg de key-operator om het sys- teem te spoelen en het systeem op lek- kage te controleren.
	Puntkegelafdichtingen of pakkin- gen oud of defect.	Raadpleeg de paragraaf <b>"Klantsupport"</b> [▶ 203]. Raadpleeg de key-operator om het sys- teem te spoelen en het systeem op lek- kage te controleren.
	Pipetteerkop is defect.	Raadpleeg de paragraaf <b>"Klantsupport</b> " [▶ 203].
Enkel kanaal lekt	Puntkegelafdichting, pakkingen of andere afdichtingen in de pipet- teerkop zijn defect.	Raadpleeg de key-operator om het sys- teem te spoelen en het systeem op lek- kage te controleren. Raadpleeg de paragraaf <b>"Klantsupport"</b> [▶ 203].
Enkele wegwerppunt niet correct opgeno- men	Individuele wegwerppunt is defect.	Vervang wegwerppunten.
	Puntkegelafdichting op deze weg- werppunt-positie is defect.	Probleem kan niet worden verholpen. Raadpleeg de paragraaf <b>"Klantsupport"</b> [▶ 203].
Enkele wegwerppunt niet gevallen	Individuele wegwerppunt is defect. Puntkegelafdichting op deze weg- werppunt-positie is defect.	Probleem kan niet worden verholpen. Raadpleeg de paragraaf "Klantsupport" [> 203].
	Onjuiste vochtigheid	Zorg ervoor dat de vochtigheid binnen de vochtigheidsgrenzen voor gebruik is. Zie paragraaf "Omgevingsvoorwaarden" [> 45].
Meerdere of alle weg- werppunten niet ge- vallen	Verkeerde wegwerppunten ge- bruikt.	Gebruik altijd wegwerppunten van Te- can. Probleem kan niet worden verholpen. Raadpleeg de paragraaf "Klantsupport"



Probleem/fout	Mogelijke oorzaak	Correctiemaatregel
Wegwerppunt-box beweegt met weg- werppunten omhoog bij het opnemen van wegwerppunten	Drager is niet correct afgesteld. X en/of Y-offsets onjuist gespeci- ficeerd.	<ul> <li>Stel alle dragers (mechanisch) precies af.</li> <li>Vervang de wegwerppunt-drager.</li> <li>Probleem kan niet worden verholpen.</li> <li>Raadpleeg de paragraaf "Klantsupport"</li> <li>[▶ 203].</li> </ul>
	Wegwerppunt-box voldoet niet aan specificaties.	Gebruik altijd wegwerppunt-boxen die voldoen aan de standaards van de "So- ciety of Biomolecular Screening". Probleem kan niet worden verholpen. Raadpleeg de paragraaf "Klantsupport" [▶ 203].
	Wegwerppunt-drager is defect (storing van wegwerppunt-box- houders).	Gebruik altijd wegwerppunt-boxen die voldoen aan de standaards van de "So- ciety of Biomolecular Screening". Probleem kan niet worden verholpen. Raadpleeg de paragraaf "Klantsupport" [▶ 203].
Onnauwkeurige pipet- teerresultaten	Wegwerppunten worden niet cor- rect opgenomen. Vloeistofverwerkingsparameters zijn niet correct. Dragers zijn niet correct afgesteld. Pipetteerkop is defect.	Raadpleeg de key-operator om het toe- passingsscript en de dragers te contro- leren. Raadpleeg de key-operator om de om- gevingsparameters en de afgiftehoogte te controleren. Controleer de puntcoating. Indien nodig, het puntblok vervangen.
	Coating van standaardpunten is beschadigd.	Raadpleeg de key-operator om het toe- passingsscript en de dragers te contro- leren. Controleer de puntcoating. Indien nodig, het puntblok vervangen.
Residu-overdracht	Coating van standaardpunten is beschadigd.	Controleer coating. Vervang vastepunt- blok of vastepuntadapter.
	Ongeschikt toepassingsscript.	Raadpleeg de key-operator om de was- of vloeistofverwerkingsinstellingen te controleren en een andere wasbuffer te gebruiken.
	Waskanalen zijn verstopt.	Reinig wasstation. Probleem kan niet worden verholpen. Raadpleeg de paragraaf "Klantsupport" [> 203].



## 8.2.5 Robotgrijperarm (RGA) verhelpen van storingen

Tab. 32: Tabel voor het verhelpen van storingen robotgrijperarm

Probleem/fout	Mogelijke oorzaak	Correctiemaatregel
Multititerplaat niet ge- pakt	Geen multititerplaten op drager. Grijpervingers kunnen de multiti- terplaat niet opnemen.	Plaats multiterplaat op drager. Stel grijperpositie in. Reinig RGA-grijpervingers.
Ongebruikelijk geluid tijdens armbeweging	Onderdelen beschadigd of versle- ten.	Raadpleeg de paragraaf "Klantsupport" [▶ 203].
Excentrische grijper- vingers onjuist uitge- lijnd	Reservevingers botsen. Vingerschroeven niet vast genoeg aangedraaid.	Lijn de excentrische grijpervingers uit. Zie paragraaf "Grijpervingeruitlijning con- troleren" [> 191].
		Gebruik een momentschroevendraaier om de schroeven aan te draaien met 3 Nm, zoals beschreven in paragraaf <b>"Ba- sis grijpervingeruitlijning voor FES-grijper-</b> <b>vingers"</b> [▶ 192].

#### 8.2.5.1 Robotgrijperarm met lange Z-as (RGA-Z) verhelpen van storingen

Tab. 33: Tabel voor het verhelpen van storingen van robotgrijperarm met lange Z-as

Probleem/fout	Mogelijke oorzaak	Correctiemaatregel
Multititerplaat niet ge- pakt	Geen multititerplaten op drager. Grijpervingers kunnen de multiti- terplaat niet opnemen.	Plaats multiterplaat op drager. Stel grijperpositie in. Reinig RGA-grijpervingers.
	Grijpervingers zijn glad.	Reinig RGA-grijpervingers.
Ongebruikelijk geluid tijdens armbeweging	Onderdelen beschadigd of versle- ten.	Raadpleeg de paragraaf "Klantsupport" [> 203].

### 8.2.6 Verhelpen van storingen in wassysteem

Tab. 34: Tabel voor het verhelpen van storingen in wassysteem

Probleem/fout	Mogelijke oorzaak	Correctiemaatregel
Foutieve overstroom- en leegloopfouten tij- dens procescyclus	Overstroomsensor niet schoon. Sensors zijn niet aangesloten of zijn defect.	Controleer of de sensor is aangesloten. Zie de handleiding van de sensorfabri- kant.
Regelmatig overstro- men van wasstation	Afvoerpomp is defect. Afvoerpompbuizen zijn geknikt, verstopt of beschadigd.	Controleer afvoerbuis. Vervang afvoerbuis indien nodig. Raad- pleeg de paragraaf "Klantsupport" [> 203].



Probleem/fout	Mogelijke oorzaak	Correctiemaatregel
Geen wasvloeistof door het wasstation gepompt of het was- systeem is leeg	Afvoerpompbuizen zijn geknikt, verstopt of beschadigd.	Controleer afvoerbuis. Vervang de afvoerpomp indien nodig. Raadpleeg de paragraaf "Klantsupport" [> 203].
	Wasstation is niet aangesloten. Wasreservoir(s) leeg of ontbreken.	Vul of vervang wasreservoir(s). Sluit het wasstation (MCA) correct aan. Zie paragraaf "Wasstation aansluiten (MCA 384)" [> 170].
	Afvoerpomp is defect.	Controleer afvoerpomp. Vervang de afvoerpomp indien nodig. Raadpleeg de paragraaf "Klantsupport" [> 203].
Wasstation over- stroom	Wasbuis is onder vloeistofopper- vlak in het afvoerreservoir.	Gebruik een wasreservoir met vaste wasbuisinlaat.
	Wegwerppunten of algen blokke- ren het wasstation.	Reinig het wasstation. Zie paragraaf "Het systeemvloeistofreser- voir en afvoerreservoir reinigen" [> 146].
	Afvoerbuis is geknikt.	Controleer buizen op knikken. Zie paragraaf "Buizen van systeemvloei- stofreservoir en afvoerreservoir controle- ren" [> 93].



## 8.2.7 Fluent ID verhelpen van storingen

Tab. 35: Tabel voor het verhelpen van storingen van Fluent ID

Probleem/fout	Mogelijke oorzaak	Correctiemaatregel
Barcode niet gelezen	Barcodelabel niet tegenover de scanner.	Laad de buisrunner uit, draai de buizen zodanig dat de barcodelabel naar links gericht zijn. Herlaad de buisrunner van de Fluent.
	Runner te snel ingeladen.	Laad de buisrunner uit en laad hem op- nieuw in.
	Slechte labelkwaliteit.	Voer de barcode handmatig in, of infor- meer de key-operator over het pro- bleem.
	Scannervenster is vuil.	Reinig het scannervenster. Zie paragraaf "Wekelijkse systeemverzorging" [> 123].
	Reflector is vuil.	Reinig de reflector. Zie paragraaf "We- kelijkse systeemverzorging" [▶ 123].
	Barcodetype of barcodelengte voor de methode niet vooraf ge- definieerd.	Informeer de key-operator over het pro- bleem.
Buisaanwezigheid niet gedetecteerd	Barcodelabelpositie te laag op de buis.	Informeer de key-operator over het pro- bleem.

## 8.2.8 Software verhelpen van storingen

Tab. 36: Tabel voor het verhelpen van storingen van software

Probleem/fout	Mogelijke oorzaak	Correctiemaatregel
Gebruikeraanmeld- scherm niet weerge- geven als dit wordt verwacht.	Gebruikersbeheer is niet geacti- veerd in FluentControl.	Raadpleeg de key-operator om het ge- bruikersbeheer te activeren.
Gebruiker kan zich niet aanmelden.	Wachtwoord is onjuist of account is geblokkeerd.	Raadpleeg de key-operator om het wachtwoord of de account te resetten.
Niet alle serviceacties voltooid. Waarschu- wing verschijnt elke keer, als FluentCon- trol wordt gestart.	Niet alle beoogde serviceacties zijn geselecteerd als voltooid in de instrumentconfiguratie.	Raadpleeg de paragraaf <b>"Klantsupport"</b> [▶ 203].



Probleem/fout	Mogelijke oorzaak	Correctiemaatregel
Touchscreen reageert niet op aanraken.	Software-stuurprogramma niet geïnstalleerd.	Raadpleeg de computerbeheerder over installatie van de stuurprogramma's op de installatie-cd en configuratie van het touchscreen.
	Touchscreen-interface is niet cor- rect geconfigureerd.	Open de instellingen van het stuurpro- gramma van het touchscreen en zorg ervoor dat het touchscreen correct is gemapt.
Touch-interface wordt niet weergegeven op het touchscreen.	Touchscreen was niet aan, toen de software werd gestart.	Schakel het instrument in en herstart de software of controleer de Touch Tool- instellingen in het FluentControl-confi- guratiesysteem.
Fout bij het starten van FluentControl.	FluentControl (SystemSW.exe) is al actief op de achtergrond (Taakbeheer).	Open Taakbeheer, bewerk Sys- temSW.exe en herstart FluentControl. Of herstart de computer.
FluentControl com- municeert niet met aangesloten hardw- areapparaten.	FluentControl is niet correct ge- configureerd voor communicatie met hardwareapparaten.	Raadpleeg de voor systeemconfiguratie verantwoordelijke persoon om de I/O- status van de hardwareapparaten acti- veren.





## 8.3 Verhelpen van storingen handelingen

## 8.3.1 Wasstation aansluiten (MCA 384)

Sluit het MCA 384-wasstation van het wassysteem als volgt aan:

- 1. Schakel het instrument uit.
- 2. Sluit de wasbuis (A) en de afvoerbuis (B) aan op het wasstation.



3. Sluit de wasbuis (A) en de afvoerbuis (B) aan.







4. Sluit alle buizen (C) aan op de wasbesturingseenheid van de MCA 384.

5. Raadpleeg de key-operator over de bedrijfsgereedheid.

## 8.3.2 Laboratoriummateriaal positioneren

Positioneer laboratoriummateriaal voor exacte armtoegang als volgt in het nest:

1. Plaats het laboratoriummateriaal (A) op het nest (B).







2. Duw het laboratoriummateriaal voorzichtig tegen de statische positioneerinrichting (C).

- 3. Schuif de diagonale schuifpositioneerinrichting (E) naar of van het laboratoriummateriaal om dit precies te laten passen.
- 4. Schuif de verticale schuifpositioneerinrichtingen (D, F) naar of van het laboratoriummateriaal om dit precies te laten passen.
- 5. Til het laboratoriummateriaal op van het nest.

Zorg ervoor dat er geen frictie is bij het plaatsen of verwijderen van de plaat.

#### 8.3.3 Inlinefilter controleren (Air FCA)

Op elk kanaal is een besturingssysteem geïnstalleerd om de Air FCApipetteerkanalen te beschermen tegen overmatig opzuigen van vloeistof.

- ✓ Het Air FCA-routineonderhoud uit methode omvat een inlinefiltercontrole die natte, beschadigde, onjuist gepositioneerde en ontbrekende filters detecteert.
- 1. Voer de **Air FCA-routineonderhoud uit** methode uit om het inlinefilter in de DiTi-kegel van een Air FCA-pipetteerkanaal te controleren.

Bij een fout moet het inlinefilter worden vervangen. Zie paragraaf "Inlinefilter vervangen (Air FCA)" [▶ 172].

#### 8.3.4 Inlinefilter vervangen (Air FCA)

Vervang het inlinefilter als volgt:

- ✓ Verontreinigde wegwerppunt-kegel.
- ✓ DiTi-kegel uit het kanaal verwijderd. Zie paragraaf "DiTi-kegel verwijderen (Air FCA)" [▶ 174].





- 1. Doorprik het inlinefilter (B) aan de zijkant met filter-verwijderingsgereedschap (C).
- 2. Wrik het inlinefilter eruit met het filter-verwijderingsgereedschap. Het filter kan zijn verontreinigd met procesvloeistoffen.
- 3. Voer het inlinefilter af.



4. Reinig de DiTi-kegel (A) met alcohol.

DiTi-kegel moet droog zijn voor hermontage.

- 5. Plaats het nieuwe inlinefilter op een schoon en effen oppervlak.
- 6. Duw het inlinefilter in de DiTi-kegel. Het inlinefilter mag niet uit de DiTi-kegel steken.
- 7. Controleer het inlinefilter volgens de door uw key-operator gedefinieerde methode.



## 8.3.5 DiTi-kegel verwijderen (Air FCA)

Verwijder de DiTi-kegel (Air FCA) als volgt:

- ✓ DiTi-steeksleutel beschikbaar.
- 1. Schakel het instrument uit.
- 2. Open het veiligheidspaneel aan de voorzijde.
- 3. Til alle Z-stangen in hun hoogste positie.
- 4. Beweeg alle Z-stangen richting de voorzijde van het instrument.
- 5. Spreid de Z-stangen zo breed als mogelijk.
- 6. Houd de puntadapter (D) en de punt-uitwerpbuis (C) vast.
- 7. Schroef de DiTi-kegel (A) los met de DiTi-steeksleutel (B).



8. Trek de DiTi-kegel voorzichtig omlaag.





In sommige gevallen is de punt-uitwerpbuis (C) of de adaptercilinder (B) nog aan de DiTi-kegel (A) bevestigd. Zie paragraaf "DiTi-uitwerpbuis monteren (Air FCA)" [> 175].



## 8.3.6 DiTi-uitwerpbuis monteren (Air FCA)

Monteer de DiTi-uitwerpbuis (Air FCA) als volgt:

- ✓ De DiTi-uitwerpbuis werd volgens de instructies verwijderd.
- ✓ DiTi-steeksleutel beschikbaar.
- 1. Steek de afdichthuls (B) in de adaptercilinder (A).





2. Schroef de DiTi-kegel (C) op de gemonteerde cilinder. Zorg ervoor dat de zwarte O-ring niet zichtbaar is, zoals weergegeven in de onderstaande afbeelding.



3. Plaats de naalduitstoterbuis (D) met de niet-kerfzijde in de gemonteerde cilinder zoals hieronder afgebeeld.



## 8.3.7 DiTi-kegel installeren (Air FCA)

Installeer de Air FCA-DiTi-kegel als volgt:

- ✓ DiTi-kegel is succesvol gemonteerd: Zie paragraaf "DiTi-uitwerpbuis monteren (Air FCA)" [▶ 175].
- ✓ DiTi-steeksleutel beschikbaar.
- 1. Schuif de adaptercilinder in de punt-uitwerpbuis (C).
- 2. Houd de puntadapter (D) en de punt-uitwerpbuis (C) vast.





3. Schroef de DiTi-kegel (A) erin met de DiTi-steeksleutel (B).

4. Voer de Air FCA-routineonderhoud uit methode.



## 8.3.8 De DiTi-optie verwijderen (FCA)

Verwijder de DiTi-optie als volgt:

- ✓ DiTi-steeksleutel
- 1. Schakel het instrument uit.
- 2. Open het frontveiligheidspaneel.
- 3. Til alle Z-stangen in hun hoogste positie.
- 4. Beweeg alle Z-stangen richting de voorzijde van het instrument.
- 5. Spreid de Z-stangen zo breed als mogelijk.
- 6. Houd de puntadapter (D) en de punt-uitwerpbuis (C) vast.
- 7. Schroef de DiTi-kegel (A) los met de DiTi-steeksleutel (B).



8. Trek de DiTi-kegel voorzichtig omlaag.

## 8.3.9 De DiTi-optie installeren(FCA)

Installeer de DiTi-optie als volgt:

- 1. Schroef de adaptercilinder (B) in de punt-uitwerpbuis (C).
- 2. Schroef de DiTi-kegel (A) op de adaptercilinder. Gebruik de DiTi-steeksleutel.







3. Duw de buis over de plastic naald tot de buis stevig op de DiTi-optie is bevestigd.

- 4. Houd de puntadapter (D) en de punt-uitwerpbuis (C) vast.
- 5. Schroef de DiTi-kegel (A) erin met de DiTi-steeksleutel (B).



## 8.3.10 Vaste punten controleren

Controleer de vaste punten:

## LET OP

## Pipetteer-onnnauwkeurigheid en fouten bij de vloeiststofdetectie!

Verbogen of beschadigde puntcoating leidt tot pipetteer-onnnauwkeurigheid en fouten bij de vloeiststofdetectie.

• Werk in geen geval met beschadigde of verbogen punten.



- 1. Schakel het instrument uit.
- 2. Open het veiligheidspaneel aan de voorzijde.
- 3. Controleer de vaste punten.
- 4. Controleer de vastepuntcoating met een spiegel.

Zorg ervoor dat vaste punten niet verbogen zijn. Als de vastepuntcoating is beschadigd of als de vaste punt is verbogen, moet deze worden vervangen. Zie paragraaf "Vaste punten verwijderen" [> 180].

#### 8.3.11 Vaste punten verwijderen

Verwijder de vaste punten als volgt:

- ✓ Vaste punten zijn gereinigd. Zie paragraaf "Systeemverzorgingstabellen"
   [▶ 117].
- ✓ Vaste punten werden gecontroleerd. Zie paragraaf "Vaste punten controleren" [▶ 179].



- E Pipetteerbuis
- 1. Schakel het instrument uit.
- 2. Open het veiligheidspaneel aan de voorzijde.
- 3. Til alle Z-stangen (A) handmatig in hun bovenste positie.
- 4. Spreid de Z-stangen zo breed als mogelijk.
- 5. Als er een afstelbare vaste punt is geïnstalleerd, de vier puntafstelschroeven losdraaien.
- 6. Schroef gelijktijdig borgmoer (C) los en houd de vaste punt met de andere hand onmiddellijk onder borgmoer vast.
- 7. Verwijder de borgmoer (C) door deze over de puntas te schuiven. *Voorkom contact tussen de borgmoer en de puntcoating.*


- 8. Als de punt (D) afstelbaar is, de borgmoer (C) omdraaien boven een schoon oppervlak om de O-ring en de sluitring te verwijderen.
- Als het kanaal is uitgerust met de klein volume-optie, de flens op de magneetklep losdraaien om de pipetteerbuis (E) die door de Z-stang (A) verloopt los te maken.
- 10. Trek de pipetteerbuis (E) een stuk (a) uit de puntadapter (B) door trekken aan de punt (D).

Gebruik een droog stuk schuurpapier voor een betere grip op de pipetteerbuisniet op de punt.

#### 8.3.12 Vaste punten installeren



Installeer vaste punten als volgt:

1. Trek pipetteerbuis voorzichtig ca. 25 mm (1 inch) (a) uit de puntadapter.

Gebruik een klein stukje schuurpapier voor een betere grip om de slang aan het eind vast te pakken.

Als een punt al eerder werd gemonteerd, met een scherp mes voor een gladde snede ca. 5 mm (0,2 inch) (b) van de buis afsnijden.





2. Bij Te-PS-punten of laag-volumepunten:

Gebruik de Te-PS-buisverbreder (G) om het buiseinde te verbreden door de Te-PS-buisverbreder tot het handvat in de buis te duwen en hierbij het gereedschap te draaien.

Als de buis nog wijd is, de Te-PS-punt ca. 4 mm (0,16 inch) in het buiseinde duwen.

3. Schuif de borgmoer op de punt.

Als de punt verstelbaar is (bijv. Te-PS), de borgmoer over de sluitring (H) en de O-ring (I) schuiven.

# *LET OP! Vermijd contact met het kwetsbare uiteinde van de punt en de coating.*

- 4. Steek de punt en de pipetteerbuis in de puntadapter.
- 5. Schroef de borgmoer op de puntadapter en draai deze vast.

Als de punt kan worden afgesteld (bijv. Te-PS), de borgmoer aandraaien tot de vier punt-afstelschroeven (K) zich in een hoek van 45° ten opzichte van het X/ Y-coördinatensysteem van het dek bevinden.

- 6. Reinig de vaste punten. Zie paragraaf "Einde van de dag" [▶ 119].
- 7. Voer een pipetteerprecisietest uit zoals gedefinieerd door de key-operator.





#### 8.3.13 Doorprikpunten verwijderen

Verwijder de doorprikpunt als volgt:

- ✓ Het instrument is uitgeschakeld.
- 1. Open het frontveiligheidspaneel.
- 2. Til alle Z-stangen in hun hoogste positie.
- 3. Beweeg alle Z-stangen richting de voorzijde van het instrument.
- 4. Spreid de Z-stangen zo breed als mogelijk.





5. Bescherm de doorprikpunten met doorprikpuntbeschermingen. Begin met de achterste doorprikpunt.

6. Schroef de borgmoer los, en houd hierbij gelijktijdig de doorprikpunt direct onder de borgmoer met de andere hand vast.





7. Trek de pipetteerbuis ca. 25 mm uit de puntadapter door aan de punt te trekken. Houd de doorprikpunt bij het trekken vast aan het bovenste uiteinde.

- 8. Trek de doorprikpunt van de buis, houd hierbij de buis met de andere hand tegen.
- 9. Verwijder de doorprikpuntbescherming niet. Voer deze samen met de doorprikpunt af in het biologische afvoerreservoir.







#### 8.3.14 Doorprikpunten installeren

- ✓ Het instrument is uitgeschakeld.
- ✓ Er is een key-operator beschikbaar.
- 1. Open het frontveiligheidspaneel.
- 2. Til alle Z-stangen in hun hoogste positie.
- 3. Beweeg alle Z-stangen richting de voorzijde van het instrument.
- 4. Spreid de Z-stangen zo breed als mogelijk.





5. Open de verpakking van de doorprikpunt. Verwijder de puntbescherming (F) niet.



Installatievolgorde van de doorprikpunten: van achter naar voren

- 6. Trek de pipetteerbuis voorzichtig ca. 25 mm uit de puntadapter.
- 7. Duw het blanke, conische einde van de doorprikpunt in het buiseinde.



8. Steek de doorprikpunt en de pipetteerbuis in de puntadapter.





9. Schroef de borgmoer op de puntadapter en draai deze handmatig vast.

- 10. Open de borgmoer iets. Beweeg de puntbescherming iets omlaag voor toegang tot de doorprikpuntschacht. Verwijder de puntbescherming nog niet helemaal.
- 11. Draai de doorprikpunt tot de puntopening tegenover de voorzijde van het instrument is. Houd de doorprikpunt met de ene hand in deze richting en draai de borgmoer met de andere hand aan.
- 12. Controleer of alle puntopeningen naar de voorzijde van het instrument gericht zijn.







13. Na installeren van alle doorprikpunten kunnen alle puntbeschermingen worden verwijderd. Begin met de achterste doorprikpunt.

- 14. Neem contact op met een key-operator om de teller in FluentControl te resetten.
- 15. Raadpleeg een key-operator voor het uitvoeren van een QC kit-controle. Zie Referentiedocumenten.
- 16. Voer de Doorprikken FCA lekkage methode.
- 17. Voer een pipetteerprecisietest (aanbeveling: QC-kit gebruiken) uit zoals gedefinieerd door de key-operator.



#### 8.3.15 Vastgeklemde doorprikpunten terugtrekken



Als een piercing tip zo vast zit dat deze niet via de software kan worden teruggetrokken, moet deze handmatig worden verwijderd.

Trek vastgeklemde piercing tips als volgt terug:

- ✓ Het instrument is uitgeschakeld.
- 1. Veiligheidspaneel aan voorzijde openen.
- 2. Til alle teruggetrokken Z-stangen handmatig in hun hoogste positie.
- 3. Bescherm alle teruggetrokken piercing tips met piercing tipbeschermingen (A). Begin met de achterste piercing tip.





4. Plaats het piercing tip-verwijderingsgereedschap naast de vastgeklemde punt op een robuust en stabiel oppervlak en breng het aan onder de borgmoer.



- 5. Draai de knop van het piercing tip-verwijderingsgereedschap tot de punt er compleet is teruggetrokken.
- 6. Draai de knop in tegengestelde richting en laat het uittrekgereedschap ongeveer 1 cm neer.
- 7. Verwijder het piercing tip-verwijderingsgereedschap. De vastgeklemde piercing tip is nu teruggetrokken.
- 8. Reinig het piercing tip-verwijderingsgereedschap met alcohol.
- 9. Controleer de piercing tip op beschadiging (bijv. verbogen piercing tip, beschadigde punt).
- 10. Vervang de piercing tip, als deze beschadigd is. Zie paragraaf "Doorprikpunten verwijderen" [▶ 183] en paragraaf "Doorprikpunten installeren" [▶ 186].
- 11. Verwijder alle piercing tipbeschermingen door het vasthouden van de borgmoer met de ene hand en verwijderen van de puntbeschermingen met de andere hand. Begin met de achterste piercing tip.
- 12. Reinig het piercing tip-verwijderingsgereedschap met alcohol.

#### Also see about this

- "Doorprikpunten verwijderen" > 183
- "Doorprikpunten installeren" > 186

#### 8.3.16 Grijpervingeruitlijning controleren

Heruitlijning van grijpervingers kan nodig zijn na een botsing of bij het implementeren van reserve-grijpervingers. Dit geldt voor alle armen in combinatie met grijpers met grijpervingers.



Onjuiste uitlijning na een botsing:

- Analyseer de situatie.
- Evalueer potentiële oorzaken van de botsing of oorzaak van de onjuiste uitlijning van de vingers, bijvoorbeeld een niet-uitgelijnde lade van een lezer, ring, onjuist geteacht/gepositioneerd hotel of een ander segment.
- Selecteer hieronder een procedure op basis van de nauwkeurigheidseisen.





- Voer een basisuitlijning uit, als er voor grijpervingers geen bovengemiddelde precisie nodig is. Zie paragraaf "Basis grijpervingeruitlijning voor FESgrijpervingers" [> 192] of paragraaf "Basis grijpervingeruitlijning voor vaste grijpervingers" [> 193].
- Als grijpervingers moeten voldoen aan specifieke eisen (Z-afwijking < ± 0,2 mm), voert u de procedure voor geavanceerde grijpervingeruitlijning uit. Zie paragraaf "Geavanceerde grijperuitlijning voor FES-grijpervingers" [▶ 194] of paragraaf "Geavanceerde grijperuitlijning voor vaste grijpervingers" [▶ 194].</li>
- 3. De grijpervingers kunnen worden gemonteerd met twee verschillende schroeven:
  - a) Torx-schroef M4x12, aangedraaid met een aanhaalmoment van 3 Nm.



b) Inbusschroef M4x12 in combinatie met een spanring (let op de positie volgens onderstaande illustratie), aangedraaid met een aanhaalmoment van 3,5 Nm.





Als er geen momentschroevendraaier beschikbaar is, de schroef aandraaien tot de ring platgedrukt is en de weerstand toeneemt. Draai dan nog 1/12 rotatie verder. Dit komt overeen met ongeveer 3,5 Nm.

#### 8.3.17 Basis grijpervingeruitlijning voor FES-grijpervingers

Dit geldt voor alle armen in combinatie met grijpers met grijpervingers.

Voer een basisuitlijning als volgt uit:

- ✓ Onjuiste uitlijning is duidelijk zichtbaar.
- ✓ Bovengemiddelde precisie is niet vereist.
- ✓ Momentschroevendraaier beschikbaar (met een optie voor 3 of 3,5 Nm). Als er geen momentschroevendraaier beschikbaar is: Torx-schroef: draai de schroeven stevig aan, maar zonder te forceren. Inbusschroef: zie "Grijpervingeruitlijning controleren" [▶ 191].



- 1. Verwijder grijpervinger van de grijperkop.
- 2. Maak de schroef tussen grijpervinger en FES-vingeradapter los.
- 3. Duw de grijpervinger tegen de bovenste en achterste stop van de adapter zoals afgebeeld in onderstaande afbeelding, en draai de schroef aan met een momentschroevendraaier (3 of 3,5 Nm).



#### 8.3.18 Basis grijpervingeruitlijning voor vaste grijpervingers

Voer een basisuitlijning als volgt uit:

- ✓ Onjuiste uitlijning is duidelijk zichtbaar.
- ✓ Bovengemiddelde precisie is niet vereist.
- ✓ Momentschroevendraaier beschikbaar (met een 3 of 3,5 Nm optie).
- 1. Maak de bevestigingsschroef (A) los.





2. Duw de grijpvinger tegen de bovenste en achterste stop van de grijpvingerbevestiging en draai de schroef aan met een momentschroevendraaier (3 of 3,5 Nm).

#### 8.3.19 Geavanceerde grijperuitlijning voor FES-grijpervingers

Dit geldt voor alle armen in combinatie met grijpers met grijpervingers.

- ✓ Momentschroevendraaier beschikbaar (met een optie voor 3 of 3,5 Nm). Als er geen momentschroevendraaier beschikbaar is: Torx-schroef: draai de schroeven stevig aan, maar zonder te forceren. Inbusschroef: zie "Grijpervingeruitlijning controleren" [▶ 191].
- Gebruik het bewegingsgereedschap om de Z-hoogte naar een hoogte van ca.
   3 mm boven de werktafel te bewegen.
- 2. Maak de schroef tussen grijpervinger en de FES-vingeradapter los.



3. Zorg ervoor dat de FES-vingeradapter stevig op de grijperkop is bevestigd. *De vingeradapters worden door een magneet op hun plaats gehouden.* 

- Gebruik het bewegingsgereedschap om de Z-hoogte naar een hoogte van 0 mm boven de werktafel te bewegen. Beweeg de laatste tiende van een millimeter zeer langzaam.
   Opmerking: Raadpleeg uw key-operator, als u geen toegang heeft tot het bewegingsgereedschap of als u dit niet kunt bedienen.
- 5. Duw de grijpervinger tegen de grijperkop en het referentievlak zoals afgebeeld in de afbeelding en draai de schroef aan met een aanhaalmoment van 3 of 3,5 Nm.
- 6. Controleer de afstelling door de kop met de hand naar 90°, 180° en 270° te draaien. Een onjuiste uitlijning op de verschillende posities duidt op een verkeerde uitlijning van de kop of de arm. In dit geval moet een FSE de uitlijning controleren.

#### 8.3.20 Geavanceerde grijperuitlijning voor vaste grijpervingers

✓ Momentschroevendraaier beschikbaar (met een 3 of 3,5 Nm optie).



- 1. Gebruik het bewegingsgereedschap om de Z-hoogte naar een hoogte van ca. 3 mm te bewegen.
- 2. Maak de schroef tussen grijpvinger en de grijpkop los.
- Gebruik het bewegingsgereedscap om de Z-hoogte naar een hoogte van 0 mm te bewegen.
   Beweeg de laatste tiende van een millimeter zeer langzaam.
   Opmerking: Raadpleeg uw key-operator, als u geen toegang heeft tot het
  - bewegingsgereedschap of als u dit niet kunt bedienen.
- 4. Duw de grijpvinger tegen de grijpkop en het referentievlak en draai de schroef aan met een aanhaalmoment van 3 of 3,5 Nm.
- Controleer de afstelling door handmatig roteren van de kop met 90°, 180°, 270°. Een onjuiste uitlijning van de verschillende posities duidt op een verkeerde uitlijning van de kop of de arm. In dit geval moet een technicus de uitlijning controleren.



#### 8.3.21 Segment controleren

Controleer of het segment gesloten is.



Afb. 69: Segment gesloten



Afb. 70: Segment open



#### 8.3.22 Positioneerpennen verwijderen

Verwijder de positioneerpennen als volgt:

# LET OP

#### **Botsing of procesfout!**

Botsing en andere procesfout kunnen een resultaat zijn van onnauwkeurig positioneren van elementen op een deksegment door losse positioneerpennen.

- Werk niet met de Fluent, als positioneerpennen ontbreken.
- ✓ Positioneerpennen zijn beschadigd.
- 1. Schuif de penverwijderaar over de positioneerpen.







2. Til de penverwijderaarhendel op, en trek de pen uit het deksegment.

8.3.23 Vergrendelpennen en positioneerpennen vervangen

### LET OP

#### **Botsing of procesfout!**

Botsing en andere procesfout kunnen een resultaat zijn van onnauwkeurig positioneren van elementen op een deksegment door losse positioneerpennen.

- Werk niet met de Fluent, als positioneerpennen ontbreken.
- De in de systeemverzorgingstabel aangegeven positioneerpennen zijn beschikbaar.
- 1. Verwijder segment van het dek:

Zie paragraaf "Segmenten verwijderen" [> 85].

2. Druk de nieuwe vergrendelpen (B) in de opening (A).







3. Druk de nieuwe positioneerpen (C) in de opening (A).

 Plaats segment op het dek: Zie paragraaf "Segmenten plaatsen" [▶ 84].



# 9 Verpakken, uitpakken, transport, opslag en verwijdering

Dit hoofdstuk bevat informatie over de regelgeving en verpakkingsregels die moeten worden opgevolgd.

## LET OP

Voorkom beschadiging door ongekwalificeerd en niet-bevoegd personeel!

Verpakken, uitpakken, transport en opslag mag alleen worden uitgevoerd door personeel van Tecan of personeel dat door Tecan is geautoriseerd!

• Raadpleeg de "Klantsupport" [> 203].

Voor informatie over het verplaatsen van het instrument, zie paragraaf "Het instrument op een kabinet binnen het laboratorium verplaatsen" [> 127].

# 9.1 Verpakkingslabels

Correcte en volledige markering van de verpakking helpt om onjuiste behandeling, ongevallen, verkeerde levering, gewichtsverlies en schade tijdens de opslag te voorkomen.

Symbool	Betekenis	Beschrijving
	Recyclen	Het verpakkingsmateriaal kan worden gere- cycled. Gooi het niet weg als huishoudelijk afval. Informatie over het materiaal dat voor deze verpakking wordt gebruikt staat onder het symbool.
	Deze kant om- hoog	Zorg ervoor dat de verpakking wordt ge- transporteerd en opgeslagen met de door de pijlen aangegeven zijde naar boven. Niet kantelen.
	Droog houden	Zorg ervoor dat de verpakking tijdens het transport en de opslag niet nat wordt.
	Breekbaar	Behandel de verpakking zorgvuldig. Breek- bare artikelen in de verpakking.
*	Uit de buurt van zonlicht houden	Voorkom dat de verpakking tijdens transport en opslag aan hitte wordt blootgesteld. Be- schermen tegen sterk zonlicht.

Tab. 37: Verpakkingssymbolen



Symbool	Betekenis	Beschrijving
	Niet stapelen	Verpakkingen niet stapelen. Deze verpakking is niet ontworpen om extra gewicht te dra- gen.

## 9.2 Afvoer

Dit hoofdstuk bevat informatie over de regelgeving inzake recycling die moet worden opgevolgd.

# LET OP

#### Recycling overeenkomstig de toepasbare regelgeving!

Neem de toepasbare wetgeving voor recycling in uw land in acht.

#### 9.2.1 Lokale vereisten Europese Unie

De Europese Commissie heeft de richtlijn betreffende afgedankte elektrische en elektronische apparatuur gepubliceerd (WEEE; 2012/19/EU).

Sinds augustus 2005 zijn producenten verantwoordelijk voor het terugnemen en recyclen van elektrische en elektronische uitrusting.

Markering	Toelichting
	<ul> <li>Negatieve gevolgen van niet-correcte afvalbehandeling.</li> <li>Behandel elektrische en elektronische apparatuur niet als ongesorteerd huishoudelijk afval.</li> <li>Verzamel elektrische en elektronische uitrusting gescheiden.</li> </ul>

#### 9.2.2 Lokale vereisten Volksrepubliek China

Markering voor de beperking van het gebruik van gevaarlijke stoffen in elektronische en elektrische producten

De Chinese elektronische industrienorm SJ/T11364-2014 **Markering voor de beperking van het gebruik van gevaarlijke stoffen in elektronische en elektrische producten** vereist de markering voor de beperking van het gebruik van gevaarlijke stoffen in elektronische en elektrische producten.



In overeenstemming met de eisen van SJ/T11364-2014 worden alle elektronische en elektrische Tecan-producten die in de Volksrepubliek China worden verkocht voorzien van een markering voor de beperking van het gebruik van gevaarlijke stoffen.

Markering	Toelichting
25	Deze markering geeft aan dat dit elektronische product bepaalde gevaarlijke stoffen bevat en veilig kan worden gebruikt tijdens de periode van milieuvriendelijk gebruik, maar dat het na de periode van milieuvriendelijk gebruik in het recyclingsysteem terecht moet komen.

#### 9.2.3 Andere eisen

Markering	Toelichting
Hg	<ul> <li>Deze lamp bevat kwik</li> <li>Recyclen of weggooien, zoals vereist door de van kracht zijnde lokale wetgeving.</li> </ul>



# **10 Klantsupport**

Dit hoofdstuk beschrijft welke bestanden en informatie Tecan nodig heeft om een eerste evaluatie van een probleem uit te voeren.

## 10.1 Contacten

Neem contact op met uw lokale distributeur of importeur, of een van onderstaande adressen.

Zie ook onze homepage op internet: www.tecan.com

Tab. 38: Klantsupportcontacten

Land/regio	Adres	Telefoon/fax/e-mail	
Azië	Tecan Asia Pte Ltd. 18 Boon Lay Way, #10-106 TradeHub 21 Singapore 609966 Singapore	Telefoon Fax E-mail	+65 6444 1886 +65 6444 1836 tecan@tecan.com.sg
Australië Nieuw-Zeeland Pacifische Ei- landen	Tecan Australia Pty Ltd Unit 2, 475 Blackburn Road Mount Waverly VIC 3149 Australië	Telefoon Telefoon Fax E-mail	Gratis telefoonnummer: 1300 808 403 +61 3 9647 4100 +61 3 9647 4199 helpdesk-aus@tecan.com
Oostenrijk	Tecan Sales Austria GmbH Untersbergstrasse 1a 5082 Grödig Oostenrijk	Telefoon Fax E-mail	+43 6246 8933 256 +43 6246 72770 helpdesk-at@tecan.com
België	Tecan Benelux B.V.B.A. Mechelen Campus Schaliënhoevedreef 20A 2800 Mechelen België	Telefoon Fax E-mail	+32 15 42 13 19 +32 15 42 16 12 tecan-be@tecan.com
China	Tecan (Shanghai) Laboratory Equip- ment Co., Ltd. 1F, T 15-4, #999, Ningqiao Road, Pilot Free Trade Zone, Shanghai, PRC, 201206	Telefoon E-mail	+86 40 0821 38 88 helpdesk-cn@tecan.com



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



Land/regio	Adres	Telefoon/fax/e-mail	
Spanje Portugal	Tecan Ibérica Instrumentación S.L. C/ Lepanto 151 Bajos E-08013 Barcelona Spanje	Telefoon E-mail	+34 93 595 25 31 helpdesk-sp@tecan.com
Zwitserland	Tecan Schweiz AG Seestrasse 103 8708 Männedorf Zwitserland	Telefoon Fax E-mail	+41 44 922 82 82 +41 44 922 89 23 helpdesk-ch@tecan.com
Verenigd Ko- ninkrijk	Tecan UK Ltd. Theale Court 11-13 High Street Theale, Reading, RG7 5AH Verenigd Koninkrijk	Telefoon Fax E-mail	+44 118 930 0300 +44 118 930 5671 helpdesk-uk@tecan.com
VS	Tecan US, Inc. 9401 Globe Center Drive, Suite 140, Morrisville, NC 27560 VS	Telefoon Fax Telefoon E-mail	+1 919 361 5200 +1 919 361 5201 Gratis telefoonnr. in de VS: +1 800 TECAN US of +1 800 832 2687 helpdesk-us@tecan.com
VS (Tecan Sys- tems)	Tecan Systems, Inc. 2450 Zanker Road San Jose, CA 95131 VS	Telefoon Fax E-mail	+1 408 953 3100 Gratis telefoonnummer: +1 800 231 0711 +1 408 953 3101 helpdesk-sy@tecan.com



# Afkortingen

#### ADT

Air displacement technology (air-displacementtechnologie)

#### Air FCA

Flexibele kanaalarm met luchtsysteem

#### ASM

Application Software Manual (handleiding voor toepassingssoftware)

#### CE

Conformité Européenne

#### cLLD

Capacitive Liquid Level Detection (capacitieve vloeistofniveaudetectie)

#### CNS

Common Notification System

#### DiTi

Wegwerptip

#### EMC

Electromagnetic Compatibility (elektromagnetische compatibiliteit)

#### EN

Europese norm

#### FCA

Flexibele kanaalarm

#### FES

Vingerwisselsysteem

#### GLP

Good Laboratory Practice

#### HEPA

High-Efficiency Particulate Arrestance

#### IEC

International Electrotechnical Commission

#### IQ

Installation Qualification (installatiekwalificatie)

#### ISO

International Organization for Standardization

#### LED

Light Emitting Diode

#### Liquid FCA

Flexibele kanaalarm met vloeistofsysteem

#### MCA

Multikanaalarm

#### MCA

Multikanaalarm

#### MCH

Multiple Channel Head (multikanaalkop)

#### MET

Geregistreerd handelsmerk van Eurofins EE als een nationaal erkend testlaboratorium

#### MIO

Bewaakte incubator-optie

#### MP

Microtiterplaat

# **•**TECAN•

#### NRTL

Nationaal erkend testlaboratorium

#### OM

Operating Manual (bedieningshandleiding)

#### OQ

Operating Qualification (gebruikskwalificatie)

#### Pc

Personal Computer

#### PP

Polypropyleen

#### rcf

relatieve centrifugale kracht

#### RF

Radiofrequentie

#### RGA

Robotgrijperarm

#### **RGA** lang Z

Robotgrijperarm lang hoogte

#### **RGA standard Z**

Robotgrijperarm standaard hoogte

#### RUO

Research use only (alleen voor onderzoek)

#### RWP

RapidWash Pump (snelwaspomp)

#### SN

Serienummer

#### Technicus

technicus

#### Te-Shake

Tecan-shaker

#### Te-VacS

Tecan-vacuümscheidingsring

#### USB

Universele seriële bus

#### WEEE

Waste Electrical and Electronic Equipment (afgedankte elektrische en elektronische apparatuur)

#### WRC

Was- en navulcentrum