thermo scientific



Gebruikershandleiding voor HyPerforma-mixer voor eenmalig gebruik (S.U.M.) met touchscreenconsole

DOC0042NL • Revisie C maart 2021



Inhoud

	Waarschuwingen, veiligheid en garantie-informatie	1
	Gebruik van deze handleiding	7
Hoofdstuk 1	Overzicht van de S.U.M.	9
	1.1 Inleiding	10
	1.2 Hardware-kenmerken	12
	1.2.1 Onderdelen van S.U.Mhardware	12
	1.2.2 Kenmerken van S.U.Msysteem	13
	1.3 Kenmerken van de touchscreenconsole	14
	1.3.1 Hardwareonderdelen van de touchscreenconsole	14
	1.3.2 Functies van de touchscreenconsole	14
	1.4 Controllers van derden	15
	1.5 Kenmerken van de BPC	15
	1.5.1 BPC-functie S.U.M.	15
	1.5.2 Bedrijfsdruk	16
	1.5.3 Werkvolume	16
	1.5.4 Aftappen	16
	1.5.5 Aseptische verbindingen	16
	1.5.6 Monstername	17
	1.6 Aanvullende/optionele systeemonderdelen	17
	1.6.1 Sonde-integratie	17
	1.6.2 Verplichte en optionele accessoires	18
Hoofdstuk 2	Hardwaremontage en -configuratie	22
	2.1 Werkplekvoorbereiding	23
	2.1.1 Elektrische aansluitingen	23
	2.1.2 Voorbereiding van de hardware	23
	2.2 Hardwaremontage	23
	2.2.1 Hardware uit de kist halen	23
	2.2.2 Motoreenheid voor S.U.M. van 2.000 l	24
	2.2.3 Voorbereiding van laadcellen	26
	2.2.4 Extra items voor montage	27
	2.3 Hardwareconfiguratie	30
Hoofdstuk 3	Configuratie van de touchscreenconsole	32
	3.1 De hardware van de touchscreenconsole configureren	33

	3.2 Het startscherm van de touchscreenconsole configureren	38
	3.2.1 Kennismaking met de touchscreenconsole	38
	3.2.2 Modules toevoegen aan het startscherm	39
	3.2.3 Een beheerdersprofiel maken en gebruiken	41
	3.3 Moduleoverzicht	44
	3.4 Vereiste externe onderdelen voor modules	47
	3.5 Modules configureren	47
	3.5.1 Interlocks en alarmen instellen	48
	3.5.2 De module Agitation (Roeren) configureren	48
	3.5.3 De module Mass (Massa) configureren	52
	3.5.4 De module Timer configureren	57
	3.5.5 De module BPC Pressure (BPC-druk) configureren	57
	3.5.6 De module Pumps (Pompen) configureren	60
	3.5.7 De module Fill (Vullen) configureren	62
	3.5.8 De module Harvest (Aftappen) configureren	64
	3.5.9 De pH-module configureren	67
	3.5.10 De module Conductivity (Geleidbaarheid) configureren	72
	3.5.11 De module Temperature (Temperatuur) configureren	77
	3.5.12 De module Liquid Pressure (Vloeistofdruk) configureren	84
	3.5.13 De module Auxiliary Output (Hulpuitgang) configureren	87
	3.5.14 De module Auxiliary Input (Hulpingang) configureren	89
Hoofdstuk 4	Kalibratieprocedures	93
	4.1 Pompen kalibreren	94
	4.2 pH kalibreren	95
	4.3 Geleidbaarheid kalibreren	97
	4.4 Roersnelheid kalibreren	98
	4.5 Massa kalibreren	99
	4.6 BPC-luchtdruk kalibreren	100
	4.7 Vloeistofdruk kalibreren	100
	4.8 Temperatuur kalibreren	101
Hoofdstuk 5	BPC laden en sonde inbrengen	102
	5.1 Richtliinen voor algemeen gebruik	103
	5.1.1 BPC voorbereiden en configureren	103
	5.1.2 Hanteringsinstructies BPC	103
	5.1.3 Werkvolume	103
	5.1.4 Vloeistofoverdracht	103
	5.1.5 Toediening	104

	5.2 De BPC laden	104
	5.3 De recirculatielijn configureren	109
	5.4 De BPC met lucht vullen	111
	5.5 De aandrijfas inbrengen	112
	5.6 BPC's laden voor open-top mengen	115
	5.7 Sondeaansluitingen maken	117
	5.7.1 Kleenpak-specificaties	117
	5.7.2 Ontvangst van de uitrusting	117
	5.7.3 Installatie	118
	5.7.4 Gammastraling	119
	5.7.5 Instructies voor het autoclaveren	119
	5.7.6 De verbinding tot stand brengen	120
	5.8 Sensoren en sondes inbrengen	127
	5.8.1 Temperatuursensoren inbrengen	127
	5.8.2 pH- en geleidbaarheidssondes voorbereiden	128
	5.8.3 pH- en geleidbaarheidssondes inbrengen	129
	5.9 De S.U.M. tarreren	131
Hoofdstuk 6	Bedieningsinformatie	133
	6.1 Vullen met vloeistof	134
		101
	6.2 Roeren instellen	135
	6.2 Roeren instellen 6.3 Poeder toevoegen met de Powdertainer	135 136
	6.2 Roeren instellen6.3 Poeder toevoegen met de Powdertainer6.4 De timer gebruiken	135 136 139
	 6.2 Roeren instellen 6.3 Poeder toevoegen met de Powdertainer 6.4 De timer gebruiken 6.5 Functies voor bewaking en controle 	135 136 139 139
	 6.2 Roeren instellen 6.3 Poeder toevoegen met de Powdertainer 6.4 De timer gebruiken 6.5 Functies voor bewaking en controle 6.5.1 Mengen 	135 136 139 139 139
	 6.2 Roeren instellen 6.3 Poeder toevoegen met de Powdertainer 6.4 De timer gebruiken 6.5 Functies voor bewaking en controle 6.5.1 Mengen 6.5.2 Meer poeder of vloeistof toevoegen 	135 136 139 139 139 139 140
	 6.2 Roeren instellen 6.3 Poeder toevoegen met de Powdertainer 6.4 De timer gebruiken 6.5 Functies voor bewaking en controle 6.5.1 Mengen 6.5.2 Meer poeder of vloeistof toevoegen 6.5.3 pH en geleidbaarheid aanpassen 	135 136 139 139 139 139 140 142
	 6.2 Roeren instellen 6.3 Poeder toevoegen met de Powdertainer 6.4 De timer gebruiken 6.5 Functies voor bewaking en controle 6.5.1 Mengen 6.5.2 Meer poeder of vloeistof toevoegen 6.5.3 pH en geleidbaarheid aanpassen 6.5.4 Temperatuur bijstellen 	135 136 139 139 139 140 142 143
	 6.2 Roeren instellen 6.3 Poeder toevoegen met de Powdertainer 6.4 De timer gebruiken 6.5 Functies voor bewaking en controle 6.5.1 Mengen 6.5.2 Meer poeder of vloeistof toevoegen 6.5.3 pH en geleidbaarheid aanpassen 6.5.4 Temperatuur bijstellen 6.5.5 De druk in de BPC-hoofdruimte bewaken 	135 136 139 139 139 140 142 143 144
	 6.2 Roeren instellen 6.3 Poeder toevoegen met de Powdertainer 6.4 De timer gebruiken 6.5 Functies voor bewaking en controle 6.5.1 Mengen 6.5.2 Meer poeder of vloeistof toevoegen 6.5.3 pH en geleidbaarheid aanpassen 6.5.4 Temperatuur bijstellen 6.5.5 De druk in de BPC-hoofdruimte bewaken 6.5.6 Alarmen bewaken 	135 136 139 139 139 140 142 143 144 144
	 6.2 Roeren instellen 6.3 Poeder toevoegen met de Powdertainer 6.4 De timer gebruiken 6.5 Functies voor bewaking en controle 6.5.1 Mengen 6.5.2 Meer poeder of vloeistof toevoegen 6.5.3 pH en geleidbaarheid aanpassen 6.5.4 Temperatuur bijstellen 6.5.5 De druk in de BPC-hoofdruimte bewaken 6.5.6 Alarmen bewaken 6.5.7 Andere functies bewaken 	135 136 139 139 139 140 142 143 144 144
	 6.2 Roeren instellen 6.3 Poeder toevoegen met de Powdertainer 6.4 De timer gebruiken 6.5 Functies voor bewaking en controle 6.5.1 Mengen 6.5.2 Meer poeder of vloeistof toevoegen 6.5.3 pH en geleidbaarheid aanpassen 6.5.4 Temperatuur bijstellen 6.5.5 De druk in de BPC-hoofdruimte bewaken 6.5.6 Alarmen bewaken 6.5.7 Andere functies bewaken 6.5.8 Monstername 	135 136 139 139 139 140 142 143 144 144 145 146
	 6.2 Roeren instellen 6.3 Poeder toevoegen met de Powdertainer 6.4 De timer gebruiken 6.5 Functies voor bewaking en controle 6.5.1 Mengen 6.5.2 Meer poeder of vloeistof toevoegen 6.5.3 pH en geleidbaarheid aanpassen 6.5.4 Temperatuur bijstellen 6.5.5 De druk in de BPC-hoofdruimte bewaken 6.5.6 Alarmen bewaken 6.5.7 Andere functies bewaken 6.5.8 Monstername 6.6 Aftappen 	135 136 139 139 139 140 142 143 144 144 145 146 147
	 6.2 Roeren instellen 6.3 Poeder toevoegen met de Powdertainer 6.4 De timer gebruiken 6.5 Functies voor bewaking en controle 6.5 Functies voor bewaking en controle 6.5.1 Mengen 6.5.2 Meer poeder of vloeistof toevoegen 6.5.3 pH en geleidbaarheid aanpassen 6.5.4 Temperatuur bijstellen 6.5.5 De druk in de BPC-hoofdruimte bewaken 6.5.6 Alarmen bewaken 6.5.7 Andere functies bewaken 6.5.8 Monstername 6.6 Aftappen 6.6.1 De module Harvest (Aftappen) gebruiken 	135 136 139 139 139 140 142 143 144 144 145 146 147
	 6.2 Roeren instellen 6.3 Poeder toevoegen met de Powdertainer 6.4 De timer gebruiken 6.5 Functies voor bewaking en controle 6.5 Functies voor bewaking en controle 6.5.1 Mengen 6.5.2 Meer poeder of vloeistof toevoegen 6.5.3 pH en geleidbaarheid aanpassen 6.5.4 Temperatuur bijstellen 6.5.5 De druk in de BPC-hoofdruimte bewaken 6.5.6 Alarmen bewaken 6.5.7 Andere functies bewaken 6.5.8 Monstername 6.6 Aftappen 6.6.1 De module Harvest (Aftappen) gebruiken 6.6.2 Handmatig aftappen 	135 136 139 139 139 140 142 143 144 144 145 146 147 147 149
	 6.2 Roeren instellen 6.3 Poeder toevoegen met de Powdertainer 6.4 De timer gebruiken 6.5 Functies voor bewaking en controle 6.5 Functies voor bewaking en controle 6.5.1 Mengen 6.5.2 Meer poeder of vloeistof toevoegen 6.5.3 pH en geleidbaarheid aanpassen 6.5.4 Temperatuur bijstellen 6.5.5 De druk in de BPC-hoofdruimte bewaken 6.5.6 Alarmen bewaken 6.5.7 Andere functies bewaken 6.5.8 Monstername 6.6 Aftappen 6.6.1 De module Harvest (Aftappen) gebruiken 6.2 Handmatig aftappen 6.7 Afsluiten en wegwerpen 	135 136 139 139 139 140 142 143 144 144 145 146 147 147 149 149
	 6.2 Roeren instellen 6.3 Poeder toevoegen met de Powdertainer 6.4 De timer gebruiken 6.5 Functies voor bewaking en controle 6.5 Functies voor bewaking en controle 6.5.1 Mengen 6.5.2 Meer poeder of vloeistof toevoegen 6.5.3 pH en geleidbaarheid aanpassen 6.5.4 Temperatuur bijstellen 6.5.5 De druk in de BPC-hoofdruimte bewaken 6.5.6 Alarmen bewaken 6.5.7 Andere functies bewaken 6.5.8 Monstername 6.6 Aftappen 6.6.1 De module Harvest (Aftappen) gebruiken 6.6.2 Handmatig aftappen 6.7 Afsluiten en wegwerpen 6.7.1 De S.U.M. afsluiten 	135 136 139 139 139 140 142 143 144 144 145 146 147 147 149 149 149

	6.7.3 Afvoerinformatie	151
Hoofdstuk 7	Onderhoud en probleemoplossing	152
	7.1 Richtlijnen voor onderhoud	153
	7.1.1 Richtlijnen voor routineonderhoud	153
	7.1.2 Softwareonderhoud van touchscreenconsole	154
	7.1.3 Richtlijnen voor preventief onderhoud	154
	7.2 Problemen met de BPC oplossen	155
	7.3 Problemen met de touchscreenconsole oplossen	156
Hoofdstuk 8	Specificaties en informatie over onderdelen	159
	8.1 Hardware-kenmerken	160
	8.1.1 Ontwerpkenmerken van S.U.M.'s van 50, 100 en 200 l	160
	8.1.2 Ontwerpkenmerken van S.U.M.'s van 500 en 1.000 l	161
	8.1.3 Ontwerpkenmerken van S.U.M.'s van 2.000 l	162
	8.2 Hardware-specificaties	163
	8.3 Specificaties van de touchscreenconsole	181
	8.4 BPC-specificaties	183
	8.4.1 Standaard BPC's 50 I	183
	8.4.2 Standaard BPC's 100 I	188
	8.4.3 Standaard BPC's 200 I	193
	8.4.4 Standaard BPC's 500 I	198
	8.4.5 Standaard BPC's 1.000 I	203
	8.4.6 Standaard BPC's 2.000 I	208
	8.4.7 Standaard open-top liners en mengschroefhulzen	212
	8.4.8 Aangepaste BPC-producten	214
	8.5 Specificaties van accessoires en opties	215
	8.5.1 Laadcellen	215
	8.5.2 Powdertainer-arm	217
	8.5.3 Kabelbeheersysteem	218
	8.5.4 Diverse items	219
	8.6 Configureerbare opties	224
Hoofdstuk 9	Algemene bestelinformatie	229
	9.1 Bestelinstructies	230
	9.2 Contactgegevens voor bestellen en ondersteuning	230
	9.3 Technische ondersteuning	231

Waarschuwingen, veiligheid en garantie-informatie

Gefeliciteerd! U hebt hoogwaardige Thermo Scientific[™]-apparatuur aangeschaft. We hebben de veiligheidsinformatie toegevoegd aan deze handleiding, die gebaseerd is op onze kennis en ervaring. Het is echter belangrijk dat u samenwerkt met uw afdeling Safety Management om te zorgen dat deze apparatuur in uw veiligheidspraktijken wordt geïntegreerd. Neem de tijd om uw eigen taakveiligheidsanalyse uit te voeren om alle potentiële gevaren te identificeren en beheren.



WAARSCHUWING: Zorg dat u deze gebruikershandleiding begrijpt voordat u deze apparatuur in gebruik neemt.

De Thermo Scientific[™] HyPerforma[™]-mixer voor eenmalig gebruik (S.U.M.) is ontwikkeld om gebruikt te worden onder traditionele farmaceutische omstandigheden. Een algemeen begrip van mengsystemen en hun werking is van belang voordat het systeem de eerste keer gebruikt wordt. Zorg dat u de gebruikershandleiding voorafgaand aan het gebruik hebt gelezen en begrepen. Indien u dit niet doet, kan dit leiden tot letsel en mogelijk productverlies.



WAARSCHUWING: Gevaarlijke interne spanning.

De motor van de mixer, motorregelaar en de touchscreenconsole beschikken allemaal over elektrische componenten. Er bestaat een risico op elektrische schokken en letsel. Koppel de voeding los voordat u elektrische componenten opent. Onderhoud mag alleen worden uitgevoerd door bevoegd servicepersoneel van Thermo Fisher Scientific. Thermo Fisher Scientific raadt het gebruik van standaard lock-outprocedures aan wanneer onderhoud aan elektrische componenten wordt uitgevoerd. De hoofdonderbreker van de touchscreen-console kan vergrendeld worden.



WAARSCHUWING: Er kan statische elektriciteit worden opgebouwd in BPC's.

BioProcess Containers (BPC's) kunnen als isolatie voor elektrostatische lading optreden. Indien elektrostatische lading wordt overgebracht naar een BPC kan de lading worden opgeslagen in de BPC en/of het interne product. Dit fenomeen verschilt per product en gebruikstoepassing. De eindgebruiker heeft daarom de exclusieve verantwoordelijkheid om te garanderen dat een risicobeoordeling wordt uitgevoerd en het risico op elektrostatische schokken wordt weggenomen. Waar van toepassing kan een roestvrijstalen koppeling van een productcontact geaard worden aan het frame om elektrostatische opbouw vanuit het materiaal in een BPC af te voeren. De juiste praktijk is om elektrostatische opbouw af te voeren door alle BPC's te aarden voordat ze worden aangeraakt. Wanneer met BPC's wordt gewerkt, is het gebruik van niet-geleidende materialen, zoals niet-geleidende handschoenen, aanbevolen.



WAARSCHUWING: Draaiende onderdelen-risico op verstrikking.

Draaiende en bewegende delen kunnen verwondingen veroorzaken. Houd uw handen uit de buurt van bewegende onderdelen tijdens het gebruik.

- Gebruik deze apparatuur niet tenzij de meegeleverde afscherming is aangebracht en naar behoren werkt.
- Het is de verantwoordelijkheid van de eindgebruiker om deze apparatuur te controleren en te bevestigen dat de apparatuur en veiligheidsvoorzieningen in goede staat zijn, en dat alle gebruikers getraind en zich bewust zijn van verstrikkingsgevaren en bijbehorende beschermende voorzieningen, zoals waarschuwingsborden en afschermingen.



WAARSCHUWING: Wees voorzichtig bij het gebruik van ladders en verhoogde platformen.

Voor sommige gebruikstoepassingen, zoals het laden van een BPC in een grote S.U.M., kan het gebruik van een ladder of platform vereist zijn. U dient voor te bevestigen dat de ladder gecontroleerd is en dat deze het gewicht van de gebruiker kan dragen. Wanneer u een ladder of platform gebruikt, dient u te zorgen dat deze stabiel staat, dat u drie contactpunten aanhoudt en dat de treden schoon zijn.

WAARSCHUWING: Volg de procedures voor lock-out/tag-out.

Om letsel te voorkomen, dient u de procedures voor lock-out/ tag-out van uw bedrijf te gebruiken wanneer u onderhoud aan apparatuur uitvoert, om zo elektrische, mechanische, pneumatische, hydraulische, chemische, thermische, zwaartekracht-gerelateerde en overige potentiële energie te isoleren en daardoor medewerkers te beschermen tegen het vrijkomen van gevaarlijke energie.

WAARSCHUWING: Wees voorzichtig met gevaarlijke chemische stoffen of materialen.

Personeelsleden die onderhoud uitvoeren aan apparatuur, moeten bekend zijn met de risico's van chemische stoffen of materialen die aanwezig zijn op of in de apparatuur. Maak gebruik van algemene communicatietechnieken voor risico's, zoals veiligheidsdatabladen, labels en pictogrammen om eventuele gevaren aan te geven.

WAARSCHUWING: Mogelijke besloten ruimte.

Gebruikers kunnen grotere S.U.M.-systemen betreden. Evalueer deze apparatuur aan de hand van de normen en procedures inzake besloten ruimtes.



WAARSCHUWING: Barstgevaar-lucht onder druk.

De BPC-kamer van de S.U.M. staat lichtelijk onder druk tijdens normale gebruiksomstandigheden. Normale, passieve ventilatie voorkomt dat de druk te hoog wordt in de kamer. Er moet gecontroleerd worden of de druk in de kamer en de druk in de toevoerleiding op de juiste manier zijn ingesteld.

- Inhoud onder druk
- Overschrijd een BPC-druk van 0,5 psi (0,03 bar) niet
- Overschrijd een toevoerdruk van 5 psi (0,34 bar) niet
- Controleer of het ventilatiefilter goed is geplaatst en naar behoren werkt



WAARSCHUWING: Heet oppervlak – niet aanraken.

De verwarmingsmantel is bedoeld om de binnenwand van het vat te verwarmen. Normale gebruiksomstandigheden genereren warmte en kunnen deze hete oppervlakken creëren.

- Heet intern oppervlak
- Contact met oppervlakken kan brandwonden veroorzaken
- Niet aanraken tijdens gebruik



WAARSCHUWING: Beknellingsgevaar.

De hanger van de Powdertainer op de S.U.M. kan handmatig verhoogd en verlaagd worden. U moet voorzichtig zijn om te voorkomen dat een gebruiker bekneld raakt of schade aan de apparatuur of de BPC wordt veroorzaakt.



WAARSCHUWING: Beknellingsgevaar.

Wees voorzichtig wanneer u de knijpventielen opent of sluit om te voorkomen dat een gebruiker bekneld raakt of verwond wordt.



WAARSCHUWING: De Thermo Scientific HyPerforma-mixer voor eenmalig gebruik mag niet worden geïnstalleerd in een potentieel explosieve omgeving, zoals vastgelegd in de toepasselijke EU ATEX-richtlijn. Het is de verantwoordelijkheid van de eindgebruiker om de mogelijke gevaren te beoordelen en begrijpen, zoals vermeld in de ATEX-richtlijnen 2014/34/EU.

Randaarde

Randaarde moet worden geverifieerd voordat de S.U.M. in een stopcontact wordt gestoken. Zorg dat het stopcontact goed geaard is.

Omgeving

- In bedrijf: 17 tot 27 °C, 20 tot 80% relatieve luchtvochtigheid, zonder condensvorming
- Opslag: -25 tot 65 °C
- Installatiecategorie II (overspanning) volgens IEC 664
- Hoogtelimiet: 2.000 meter

Elektrische aansluitingen

De apparatuur moet worden gevoed door een stroomkring van 15 A zonder aardlekschakelaar. Aardingsstoringen doen zich voor wanneer stroom ergens lekt en elektriciteit dus naar de grond ontsnapt. Wanneer het menselijk lichaam optreedt als het pad voor lekkage naar de grond, kan dit leiden tot elektrocutie. Een aardlekschakelaar (GFCI) controleert de stroom die naar de grond loopt en schakelt de voeding uit (activeert de GFCI) in een fractie van een seconde voor stroomspanning die veel lager is dan als gevaarlijk wordt beschouwd. Vanwege de gevoeligheid van aardlekschakelaars bij een elektrische lekkage (van een paar mA), raden we u aan de mixer voor eenmalig gebruik NIET in een stopcontact met GFCI te steken.

Informatie over de watermantel van het vat

Het S.U.M.-hardwaresysteem met watermantel is ontwikkeld om gebruikt te worden met water als hulpmiddel voor warmteoverdracht. De temperatuur mag 50 °C (122 °F) niet overschrijden bij een bedrijfsdruk van minder dan 150 psig (1 MPa). Om de veiligheid te garanderen, raden we u aan om de S.U.M. te gebruiken bij 75 psig of minder.

Opmerking: De BPC-gebruikslimieten van de S.U.M. zijn 5 tot 40 °C wat betreft temperatuur. De interne druk mag de 0,5 psi niet overschrijden. De watermantel hoeft niet geregistreerd, geïnspecteerd en van het Code U-symbool voorzien te worden volgens artikel U-1(c)2(f) van de ASME Boiler and Pressure Vessel Code en/of de Europese richtlijn 97/23/EG betreffende drukapparatuur (PED). Op aanvraag kan een conformiteitsverklaring voor goede technische PED-praktijken beschikbaar worden gesteld.

Garantie-informatie

Eventuele garantie voor deze apparatuur is niet van toepassing op: (a) normale sliitage; (b) ongelukken, rampen of overmacht; (c) verkeerd gebruik, fouten of nalatigheid door uw toedoen; (d) het gebruik van de apparatuur op een manier waarvoor het niet ontworpen is; (e) externe aantasting van de apparatuur, zoals, maar niet beperkt tot, externe doorboring, stroomuitval of elektrische piekspanningen; (f) onjuiste opslag en onjuist gebruik van de apparatuur; (g) het gebruik van de apparatuur in combinatie met apparatuur of software die niet door ons is geleverd; (h) apparatuur die aan u verkocht is als 'gebruikte' producten; (i) contact met verkeerd gebruikte of niet-goedgekeurde chemische stoffen of materialen; (j) installatie, verwijdering, gebruik, onderhoud, opslag of hantering op een ongepaste, onjuiste of nietgoedgekeurde manier, zoals, maar niet beperkt tot, het niet volgen van de meegeleverde documentatie of instructies of die gerelateerd zijn aan de apparatuur, het gebruik buiten de omgevings- of andere toepassingsspecificaties, of het gebruik in combinatie met nietgoedgekeurde software, materialen of overige producten; (k) productie in overeenstemming met de vereisten die u hebt opgegeven; (I) installatie van de software, interface of producten die niet door ons zijn goedgekeurd; (m) het gebruik van meegeleverde instructies of documentatie om wettelijke goedkeuringen te ondersteunen; (n) de prestaties, efficiëntie of compatibiliteit van vermelde componenten; en (o) de prestaties van aangepaste apparatuur, producten of vermelde componenten, of het verkrijgen van resultaten via de apparatuur, vermelde componenten of services binnen het bereik dat u had gewenst, zelfs als die bereiken aan ons zijn gecommuniceerd en beschreven in specificaties, een offerte of een arbeidsverklaring. BOVENDIEN ZULLEN BIJ ELK(E) INSTALLATIE, ONDERHOUD, REPARATIE, SERVICE, VERPLAATSING, AANPASSING OF ENIGE ANDERE VORM VAN MANIPULATIE VAN DE APPARATUUR DOOR EEN NATUURLIJKE PERSOON OF RECHTSPERSOON ANDERS DAN WIJZELF ZONDER ONZE VOORAFGAANDE SCHRIFTELIJKE GOEDKEURING, EN BIJ ELK GEBRUIK VAN VERVANGENDE ONDERDELEN DIE NIET DOOR ONS ZIJN GELEVERD, ALLE GARANTIES OP DE BETREFFENDE APPARATUUR DIRECT VERVALLEN EN WORDEN INGETROKKEN. INDIEN DE APPARATUUR IN DE VERENIGDE STATEN WORDT GEBRUIKT, KUNNEN WE UW GARANTIE INTREKKEN INDIEN U DE APPARATUUR BUITEN DE VERENIGDE STATEN VERZENDT.

Gebruiksbeperkingen

U moet deze apparatuur in overeenstemming met onze documentatie gebruiken, en, indien van toepassing, in combinatie met onze andere bijbehorende instructies, inclusief, zonder beperking, een productlabel voor 'alleen voor onderzoeksdoeleinden' of een labellicentie voor 'beperkt gebruik'. Deze apparatuur is bedoeld voor onderzoeksdoeleinden of verdere productie van bioverwerkingstoepassingen en niet voor diagnostisch gebruik of directe toediening in mensen of dieren. We leggen de apparatuur niet voor regelmatige controle voor aan een overheidsinstantie of andere organisatie en we valideren de apparatuur niet voor klinisch of diagnostisch gebruik, voor veiligheid en effectiviteit, of voor overig specifiek gebruik of toepassing.

Richtlijnen voor seismische activiteit

De koper van de apparatuur is er verantwoordelijk voor om aan de landspecifieke wetten te voldoen en voor de evaluatie van seismische waarden om te controleren of de installatie en veiligheid van de apparatuur geschikt zijn voor de betreffende locatie. De koper draagt tevens de verantwoordelijkheid voor het evalueren van de structuur van het gebouw voor de betreffende apparatuur, om te verzekeren dat de verankering en de bevestiging van de apparatuur en het gebouw bestand zijn tegen seismische activiteit. We raden de koper ten zeerste aan om een plaatselijk, bevoegd, extern architecten- en ingenieursbureau om een correcte technische analyse en gecertificeerde documentatie te vragen voorafgaand aan de installatie van de apparatuur in het gebouw. De koper is bovendien verantwoordelijk voor het vastzetten en verankeren van de apparatuur op een aangegeven, vaste locatie. Thermo Fisher Scientific kan op aanvraag en tegen overeengekomen kosten ondersteuning bieden bij de seismische verankering en bevestiging voor aangeschafte apparatuur, op basis van de bouw- en landspecifieke wetten.

Verplaatsbare apparatuur (d.w.z. niet-vaste apparatuur of op een bevestiging met zwenkwielen) komt niet in aanmerking voor seismische ontwerpvereisten in overeenstemming met ASCE 7-16, hoofdstuk 13, artikel 1.4. Hoewel deze apparaten niet in aanmerking komen voor de seismische ontwerpvereisten van ASCE 7, dient u er rekening mee te houden dat dergelijke apparatuur kan omkiepen bij seismische activiteit. Het is daarom de verantwoordelijkheid van de klant om de seismische veiligheid voor verplaatsbare apparatuur in acht te nemen op de betreffende locatie.

Gebruik van deze handleiding

Reikwijdte van deze publicatie

Deze gebruikershandleiding bevat informatie over het standaard HyPerforma S.U.M.-systeem van Thermo Scientific, met inbegrip van hardware, onderdelen, controlemethodes voor productontwerp, installatie, bediening en specificaties Ze is bestemd voor personen met of zonder ervaring met systemen van Thermo Scientific, maar die enige kennis hebben van bioproductieprocessen en grootschalige mengsystemen.

Rev.	Datum	Sectie	Aangebrachte wijziging	Auteur
А	10/2018		Oorspronkelijke versie	E. Hale
В	01/2019	8,2	'Relatieve vochtigheid' gecorrigeerd in specificaties voor alle systeemgroottes	E. Hale
В	01/2019	8,3	Sensorbereiken van laadcellen voor S.U.M.'s van 200, 500, 1.000 en 2.000 I toegevoegd aan tabel 8.19	E. Hale
В	01/2019	8,2	Algehele breedte 500 I gecorrigeerd in specificaties tot 117,4 cm (46,2 inch)	E. Hale
В	01/2019	8,2	Bereik mengsnelheden 2.000 I gecorrigeerd in specificaties tot 30–350 RPM	E. Hale
В	01/2019	8,4	Grootte in tabel 8.48 gecorrigeerd tot 2.000 I	E. Hale
В	01/2019		Kleine opmaakwijzigingen aangebracht	E. Hale
С	10/2019	8,2	Kleine revisies en lengte van wagentje bijgewerkt in afbeelding 8.8	T. Golightly

Informatie over documentwijzigingen

Vragen over deze publicatie

Indien u vragen of opmerkingen heeft over de inhoud van deze publicatie, neem dan contact op met **technicaldocumentation@ thermofisher.com** of met uw Thermo Fisher Scientific-verkoopteam.

Gerelateerde publicaties

Neem contact op met uw plaatselijke verkoopvertegenwoordiger voor informatie over de hieronder vermelde gerelateerde publicatie.

Publicatie	Beschrijving
Validatiehandleiding voor HyPerforma S.U.M. met touchscreenconsole (DOC0067)	Informatie over validatieprocedures
HyPerforma S.U.M. Gegevensbladen	Productbeschrijvingen en bestelinformatie
Integratorhandleiding touchscreen-console (DOC0069)	Informatie over de integratie van de touchscreenconsole in controllers van derden
Uitpakhandleiding voor HyPerforma S.U.M. met touchscreenconsole (DOC0061)	Instructies om het S.U.Msysteem uit te pakken

Afkortingen

Zie onderstaande lijst voor definities van afkortingen die in deze publicatie worden gebruikt.

BPC	BioProcess Container
DO	Opgeloste zuurstof
E-Stop	Noodstopknop
ETP	Specificatiepakket van de apparatuur
GFCI	Aardlekschakelaar
ID	Binnendiameter
IEC	Internationale elektrische code
OD	Buitendiameter
PED	Richtlijn drukapparatuur
P&ID	Proces- en instrumentendiagram
RTD	Weerstandtemperatuurdetector
S.U.M.	Mixer voor eenmalig gebruik
TCU	Temperatuurregelaar
VFD	Aandrijving met variabele frequentie

Overzicht van de S.U.M.

Inhoudsopgave

- 1.1 Inleiding
- 1.2 Hardware-kenmerken
- 1.3 Kenmerken van de touchscreenconsole
- 1.4 Door eindgebruiker geleverde onderdelen
- 1.5 Kenmerken van de BPC
- 1.6 Aanvullende/optionele systeemonderdelen

1.1 Inleiding

De HyPerforma mixer voor eenmalig gebruik (S.U.M.) van Thermo Scientific is een eenmalig bruikbaar alternatief voor traditioneel mengen in geroerde tanks. Het systeem is gebaseerd op hetzelfde mengprincipe als de Thermo Scientific HyPerforma bioreactor voor eenmalig gebruik (S.U.B.). Beide systemen gebruiken een mengschroef die aan een overhead mengmotor is gekoppeld via een afgedichte lagereenheid zodat de mengschroef kan draaien zonder de integriteit van het systeem in gevaar te brengen. De S.U.M. is ontworpen voor een gesloten poeder-op-vloeistof- en vloeistof-opvloeistof-systeem met contactoppervlakken voor eenmalig gebruik en open-top mengen.

Elk S.U.M.-systeem bestaat uit de volgende onderdelen:

- **1. Roestvrijstalen buitenste ondersteuningscontainer**, beschikbaar met of zonder mantelverwarmingssysteem.
- 2. Touchscreenconsole, om meerdere sensoren en functies te bewaken en in te stellen.
- **3. BioProcess Container (BPC)**, die gammabestraald wordt geleverd.



Afbeelding 1.1. S.U.M.-systeem van 100 l met alle beschikbare opties.

De **buitenste ondersteuningscontainer** is een roestvrijstalen vat dat de BPC bevat en ondersteunt. De container is ontwikkeld en gefabriceerd om elke BPC te ondersteunen en biedt gemakkelijke toegang voor bediening. De aandrijfas is verwijderbaar en herbruikbaar binnen de gestelde gebruiksparameters en wordt via de mengeenheid en in de lagerpoort aangebracht in closed-top BPC's Laadcellen zijn voor alle systemen beschikbaar om wegen te vergemakkelijken.

Met de **touchscreenconsole** kunnen verschillende functies tijdens het gebruik direct worden benut, zoals temperatuur, pH, geleidbaarheid, roersnelheid, vullen en aftappen. Dankzij het flexibele ontwerp kunnen gebruikers alarmen met toegewezen vergrendelingen inschakelen, verbinding maken met externe apparaten (zoals pompen) en deze instellen, en het startscherm aanpassen met de gewenste functionaliteit.

De **BPC** biedt gebruiksklare contactoppervlakken voor eenmalig gebruik, met inbegrip van een mengschroef, een afgedichte lagereenheid en slangen voor vloeistofoverdracht. De opties zijn:

- Open-top liner voor media-/bufferbereiding
- Gesloten BPC's met poederpoort, ontworpen voor integratie in de Thermo Scientific Powdertainer en bedoeld voor ergonomische, ingesloten media-/bufferbereiding
- Gesloten BPC's met mogelijkheden voor bewakingssondes
- Gesloten BPC's voor vloeistof-op-vloeistof mengen van kritieke steriele oplossingen

In deze gebruikershandleiding vindt u informatie over de configuratie, de bediening, het onderhoud en het oplossen van problemen voor alle S.U.M.-systemen in de volgende volumes: 50, 100, 200, 500, 1.000, en 2.000 liter.

1.2 Hardware-kenmerken

1.2.1 Onderdelen van S.U.M.-hardware

In afbeeldingen 1.2 en 1.3 hierna worden alle beschikbare onderdelen van S.U.M.-systemen in de groottes van 500 en 2.000 l geïllustreerd.



Afbeelding 1.2. Vooraanzicht van S.U.M. van 500 l.

- 1. Powdertainer-arm (optioneel)
- 2. Motor van mixer met veiligheidsklep
- 3. Gegolfde mantel 0,95 cm (3/8 inch)
- Standaard gereedschapsset: Vierkante momentsleutel van 10 mm (3/8 inch) x 16,9 Nm (150 in-lb.) en vergrendelsleutel voor laadcellen en motorkap
- 5. Buitenste ondersteuningscontainer van roestvrij staal
- 6. Aandrijfas, opgeborgen
- 7. Mand (optioneel)



Afbeelding 1.3. Vooraanzicht van S.U.M. van 2.000 I.

- 8. Touchscreenconsole
- 9. Planken (optioneel)
- 10. Wagen
- 11. Zwenkwielen (2 scharnierend, 2 vast voor S.U.M.'s van 50-1.000 l)
- 12. Haken voor kabelbeheersysteem (optioneel)
- 13. Vloeistofkijkvensters
- 14. Deur voor laden BPC
- 15. Venster sondetoegang
- 16. Sondecliphanger

1.2.2 Kenmerken van S.U.M.-systeem

De S.U.M. is ontworpen met het oog op mobiliteit, bedieningsgemak en eenvoudige integratie van disposables. **Opmerking:** De systemen van 2.000 l zijn niet draagbaar en hebben geen wielen. Hardwaretekeningen en specificatietabellen voor alle S.U.M.-groottes vindt u in hoofdstuk 8 – Specificaties en informatie over onderdelen

Schudden

De mengsnelheid van de S.U.M. kunt u instellen met de touchscreenconsole. De interface geeft de roersnelheid aan in aantal omwentelingen per minuut (RPM).

Temperatuurregulering

Het S.U.B.-hardwaresysteem met watermantel is ontwikkeld om gebruikt te worden met water als hulpmiddel voor warmteoverdracht, bij een temperatuurbereik van 2–50°C. Onder bepaalde omstandigheden kunnen hogere temperaturen mogelijk zijn. Raadpleeg uw vertegenwoordiger van Thermo Scientific als hogere temperaturen gewenst zijn. De procestemperatuur kan worden bewaakt met de temperatuurmodule van de touchscreenconsole, dankzij een temperatuursensor die wordt aangebracht in de thermowell van een BPC met sondepoorten.

Laadcellen

Laadcellen zijn beschikbaar voor alle roestvrijstalen S.U.M.-hardware. Laadcellen kunnen in de fabriek worden geïnstalleerd of later worden aangebracht door een gecertificeerde onderhoudsmonteur. Laadcellen worden ongekalibreerd geleverd. De fabrikant van de laadcellen of een gekwalificeerde monteur moet deze systemen ter plekke kalibreren.

1.3 Kenmerken van de touchscreenconsole

1.3.1 Hardwareonderdelen van de touchscreenconsole

Afbeelding 1.4 op de volgende pagina toont de hardwareonderdelen van de touchscreenconsole.



Afbeelding 1.4. Voor- en achteraanzicht van de touchscreenconsole.

1.3.2 Functies van de touchscreenconsole

Met de touchscreenconsole kunnen gebruikers meerdere functies bewaken en instellen, zoals roeren, pH, geleidbaarheid, temperatuur en massa. Op het aanpasbare startscherm kunnen de gebruikers alleen de functies weergeven die ze voor hun activiteiten nodig hebben. Zie voor meer informatie hoofdstuk 3 – Touchscreenconsole configureren

1.4 Controllers van derden

De standaard HyPerforma S.U.M.'s hebben een complete, geïntegreerde touchscreenconsole om verschillende functionaliteiten in te stellen. Controllers van derden kunnen echter worden aangesloten via de Modbus- en Profibus-poorten op de achterzijde van de touchscreenconsole. In deze configuratie wordt de touchscreenconsole gebruikt om gegevenspakketten van de controllers van derden te verzenden naar verschillende andere apparaten. Zie de Integratorhandleiding voor meer informatie. Controllers van de volgende fabrikanten zijn aanpasbaar voor gebruik met de S.U.M.:

- Thermo Fisher Scientific
- ABEC
- Bellco
- Broadley-James
- Dasgip
- Emerson
- Honeywell
- New Brunswick Scientific
- Pendotech
- Sartorius Stedim Biotech

1.5 Kenmerken van de BPC

1.5.1 BPC-functie S.U.M.

De S.U.M. BPC (een gesloten BioProcess Container of een open-top liner) bevat het mengproces. De BPC-kamer is gemaakt van CX5-14film of Aegis5-14-film. De open-top liner is gemaakt van CX3-9 film. De BPC en liner zijn gecoëxtrudeerde structuren die specifiek zijn ontworpen voor gebruik in biofarmaceutische processen. Alle materialen zijn gekwalificeerd voor een serie fysieke, mechanische, biologische en chemische compatibiliteitsvereisten. De BPC van de mixer is gammabestraald op een minimale drempelwaarde van meer dan 25 kGy. Dit resulteert in het uiteenvallen van elektronen, waardoor micro-organismen worden vernietigd of niet langer in staat zijn zich te reproduceren binnen een verpakte BPC. Het leidt echter niet tot residuen of radioactiviteit in de BPC. Er zijn twee standaard BPCconfiguraties beschikbaar voor poeder-op-vloeistof- en vloeistof-opvloeistof-toepassingen, elk verkrijgbaar met of zonder sondepoorten. De open-top liner heeft een gescheiden mengschroef en aandrijfashulssysteem voor open-top mengen Zie gedeelte 8.4 voor meer informatie over BPC's.

1.5.2 Bedrijfsdruk

VOORZICHTIG: De BPC S.U.M. is niet geclassificeerd als een drukvat. De gasdruk mag niet hoger zijn dan 0,03 bar (0,5 psi) binnen een statische BPC of 0,007 bar (0,1 psi) wanneer de motor tijdens gebruik draait,.
Opmerking: De touchscreenconsole heeft een ingeprogrammeerd instelpunt voor luchtdruk van 0,004 bar (0,06 psi). De BPC mag tijdens het opblazen niet te gespannen worden.
Overdruk kan leiden tot beschadiging van de BPC of persoonlijk letsel. De bedrijfsdruk kan worden bewaakt met de module van de BPC-druk op de touchscreenconsole.
Opmerking: Bewaking van de BPC-druk vereist een BPC met drukmonitor.

1.5.3 Werkvolume

ledere S.U.M. is ontworpen voor een specifiek werkvolumebereik. Het minimale werkvolume en het nominale werkvolume worden vermeld in de specificatietabellen in hoofdstuk 8 — Specificaties en informatie over onderdelen. De werkelijke werkvolumes mogen de aangegeven nominale werkvolumes niet overschrijden. Indien nodig kan de BPC een kleine volumeoverschrijding aan (68 I voor de S.U.M. van 50 I, 110 I voor de S.U.M. van 100 I, 220 I voor de S.U.M. van 200 I, 550 I voor de S.U.M. van 500 I, 1.100 I voor de S.U.M. van 1.000, 2.100 I voor de S.U.M. van 2.000 I). **VOORZICHTIG:** Werkvolumes kleiner dan de genoemde minima kunnen leiden tot hardwarestoringen en beschadiging van de BPC.

1.5.4 Aftappen

De S.U.M. heeft een onderste afvoerlijn waarmee vloeistof afgetapt kan worden met een peristaltische pomp of de zwaartekracht. De onderste afvoerlijn kan aangesloten worden met behulp van de meegeleverde snelaansluiting van 12,7 mm (1/2 inch). Wanneer de S.U.M. BPC wordt gehanteerd terwijl de laatste paar liter vloeistof wordt verwijderd, kan vloeistofophoping worden geminimaliseerd.

1.5.5 Aseptische verbindingen

Er zijn meerdere aseptische verbindingsopties beschikbaar voor S.U.M.-gebruikers. De standaard-BPC bevat stukken slang voor de slanglasser, snelaansluitingen voor gebruik onder een laminaire luchtstroomkap en stoombare sanitaire aansluitingen voor vloeistof-op-vloeistof-BPC's. De S.U.M. BPC is ontworpen met thermoplastische slangen van verschillende lengtes en afmetingen met het oog op toevoegen en onttrekken vanuit de S.U.M. BPC. Zie de BPC-opties in hoofdstuk 8 van deze publicatie voor aangepaste opties voor eindbehandeling.

1.5.6 Monstername

Tijdens het gebruik van de S.U.M. moeten mogelijk monsters worden genomen om verschillende, door de gebruiker ingestelde parameters te bewaken, zoals pH, spectrofotometrische analyse en osmolaliteit. Monsters kunnen op verschillende manieren uit de S.U.M. worden genomen, afhankelijk van de BPC-configuratie. Monstername is gemakkelijk uitvoerbaar via de recirculatielus en de SmartSite-poort die op alle standaard BPC-configuraties aanwezig is. In toepassingen met volledig volume waarin de recirculatielus niet wordt gebruikt, kunnen monsters direct via de poederpoort worden genomen met behulp van een van de lijnsets op de bovenkant van de BPC of via de afvoerlijn.

Voor BPC-configuraties die sondepoorten gebruiken, kan de BPC worden uitgerust met een monsterpoort voor kleine volumes die deel uit maakt van de BPC-thermowell. Met deze siliconen stijgbuis met een kleine diameter en een lengte van 15,24 cm (6 inch) kunnen monsters met een laag dood volume worden genomen. De stijgbuis wordt geleverd met een aseptische Luer-Lock-connector (SmartSite[™]) waarmee rechtstreeks monsters kunnen worden genomen of verschillende monsterverdeelstukken kunnen worden aangesloten met de standaard Luer-Lock-connector. Verdeelstukken kunnen ook met een slanglasser op de C-Flex monsterlijn worden gelast.

1.6 Aanvullende/optionele systeemonderdelen

1.6.1 Sonde-integratie

De sonde-eenheid heeft een innovatief ontwerp voor eenmalig gebruik waarin door de gebruiker geleverde pH-sondes worden verpakt en aangesloten op de S.U.M. BPC. De sonde-eenheid (afbeelding 1,5) bestaat uit de volgende onderdelen:

- 1. Gegoten balgafdekking
- 2. Sondeadapter met schroefdraad
- Pall[™] Kleenpak[™]-connector (KPCHT-serie, voor hoge temperaturen)
- 4. Kabelbinders



Afbeelding 1.5. Sonde-eenheid

Opmerking: Afbeelding 1.5 (boven) toont een sonde-eenheid met een Kleenpak aseptische connector. Uw S.U.M. BPC gebruikt in plaats daarvan mogelijk CPC[™] AseptiQuik[™] aseptische connectoren of niet-aseptische snelaansluitingen.

1.6.2 Verplichte en optionele accessoires

De volgende aanvullende accessoires zijn verkrijgbaar om het gebruik van de S.U.M. te ondersteunen. Zie hoofdstuk 8 voor meer informatie over hardwareaccessoires.

Powdertainer-hanger en -houderarm van Thermo Scientific

De Powdertainer[™]-hanger en -houderarm (afbeelding 1.6 rechts) zijn een optioneel accessoire dat wordt gebruikt om poederzakken op te hangen en te positioneren. Het is bedoeld voor S.U.M.-systemen die poeder-op-vloeistof mengen gebruiken.

Robuuste slangklemmen

Slangklemmen (afbeelding 1.7) zijn vereist om ongebruikte lijnsets handmatig dicht te knijpen en te voorkomen dat procesvloeistoffen in de lijnsets komen. Op S.U.M.-systemen met de touchscreenconsole kunnen in plaats daarvan knijpventielen worden gebruikt die automatisch worden gecontroleerd. **Opmerking:** Voordat een steriele sonde wordt ingebracht, moeten handmatige slangklemmen zijn geplaatst om sondepoorten af te sluiten. Gebruikers moeten één slangklem hebben voor elke gebruikte aansluitpoort.



Afbeelding 1.6. Robuuste slangklem.



Afbeelding 1.7. Powdertainerarm S.U.M. van 50 I.

Knijpventielen

Knijpventielen (afbeelding 1.8) worden aangesloten op de touchscreenconsole en worden gebruikt om de vloeistofstroom door de BPC-slangen automatisch te reguleren tijdens het vullen of aftappen. Het volgende bereik slangmaten kan worden gebruikt met S.U.M.-knijpventielen:

- Maximum: Slangen met een buitendiameter van 19 mm (3/4 inch) met een wand van 3,18 mm (1/8 inch)
- Minimum: Slangen met een buitendiameter van 12,7 mm (1/2 inch) met een wand van 3,18 mm (1/8 inch)



Afbeelding 1.8. Knijpventiel.

Sondeclips

Sondeclips (afbeelding 1.9) zijn vereist om sondes tijdens het gebruik van de S.U.M. op hun plek te houden.



Afbeelding 1.9. Plastic sondeclip.

Autoclaafplaat voor sondekits

De roestvrijstalen autoclaaftray (afbeelding 1.10) is ontworpen voor aseptische toepassingen en is verplicht tijdens het sterilisatieproces. De ondersteuningstray biedt een hellende bevestiging voor twee sondes tijdens het autoclaveren om de spanning op de sondes te minimaliseren en te voorkomen dat de silicone balgen worden ingedrukt. **Opmerking:** Afbeelding 1.10 laat zien hoe de autoclaaftray wordt gebruikt met Kleenpak aseptische connectoren. Voor systemen met AseptiQuik-connectoren is een andere autoclaaftray vereist. Zie paragraaf 8.5 – Specificaties van accessoires en opties voor meer informatie.



Afbeelding 1.10. Autoclaaftray.

Geleidbaarheids- en pH-sondes

Tabellen 1.1 en 1.2 tonen de lengte, de diameter het type procesverbinding, het signaaltype en de onderdeelnummers voor de geleidbaarheids- en pH-sondes die in de S.U.M. zijn geïntegreerd. De lengte- en diametervereisten zijn gebaseerd op de vereiste inbrengdiepte van de sonde bij gebruik in combinatie met de sondepoort.

Tabel 1.1. Aanbevolen geleidbaarheidssondes voor gebruik met de S.U.M.

Sonde	Onderdeelnummer Thermo Fisher	Onderdeelnummer fabrikant	Diameter	Type procesverbinding	Lengte	Type signaal
Jumo-geleiding	SV51148.01	00682581	12 mm (0,47 inch)	13,5 PG	225 mm (8,85 inch)	M12
Mettler Toledo- geleiding	SV51148.02	52001998	12 mm (0,47 inch)	13,5 PG	225 mm (8,85 inch)	VarioPin

Tabel 1.2. Aanbevolen pH-sondes voor gebruik met de S.U.M.

Sonde	Onderdeelnummer Thermo Fisher	Onderdeelnummer fabrikant	Diameter	Type procesverbinding	Lengte	Type signaal
Mettler Toledo pH	SV51147.01	59903228	12 mm (0,47 inch)	13,5 PG	225 mm (8,85 inch)	S8
Broadley James pH	SV51147.03	F-635-B225-DH	12 mm (0,47 inch)	13,5 PG	225 mm (8,85 inch)	S8
Thermo Scientific pH	SV51147.02	SV51147.02	12 mm (0,47 inch)	13,5 PG	225 mm (8,85 inch)	S8

Pompen

Pompen zijn optioneel voor S.U.M.-systemen. Indien gebruikt in combinatie met de touchscreenconsole, stellen pompen gebruikers in staat vloeistof automatisch bij te vullen en af te tappen en bolussen van zuur-, base- of andere bufferoplossingen tijdens het mengen toe te voeren. Verschillende pompmaten zijn verkrijgbaar om in te spelen op de verschillende behoeften van gebruikers.

Laadcellen

Laadcellen worden gebruikt om het gewicht van de inhoud van een S.U.M. te bepalen. Hoewel laadcellen voor alle systemen optioneel zijn, zijn ze verplicht wanneer u de modules Fill (Vullen), Harvest (Aftappen) en Mass (Massa) op de touchscreenconsole gebruikt.

Kabelbeheersysteem

Het kabelbeheersysteem is optioneel voor alle S.U.M.-systemen. Meerdere aan de buitenste ondersteuningscontainer bevestigde clips worden gebruikt om kabels van de touchscreenconsole te organiseren.

Aandrijfassen

De aandrijfas wordt via de mengeenheid en in de lagerpoort aangebracht in closed-top BPC's. In S.U.M.-eenheden van 50 en 100 I worden alleen eendelige aandrijfassen gebruikt, in eenheden van 200, 1.000 en 2.000 I alleen tweedelige. Beide typen assen zijn beschikbaar voor eenheden van 500 I.

AC- en DC-motoren

S.U.M.-systemen van 50 l, 100 l, 200 l, 500 l en 1.000 l zijn alleen verkrijgbaar met gelijkstroommotoren. S.U.M.'s van 2.000 l zijn alleen verkrijgbaar met wisselstroommotoren.

Diverse optionele items

- S.U.M.-thermowell of monsterpoort poort om de temperatuursensor te kalibreren en valideren.
- Druksensoren
- Communicatiekabels
- Sondes
- Temperatuursensor
- P&ID-tags
- Drukontlastingsklep met watermantel alleen voor S.U.M.systemen met watermantel
- Planken en mand om pompen en flessen op te zetten

Zie paragraaf 8.5 — Specificaties van accessoires en opties voor meer informatie over hardwareaccessoires.

Hardwaremontage en -configuratie

Inhoudsopgave

- 2.1 Werkplekvoorbereiding
- 2.2 Hardwaremontage
- 2.3 Hardwareconfiguratie

2.1 Werkplekvoorbereiding

2.1.1 Elektrische aansluitingen

De voedingskabel voor de touchscreenconsole kan worden besteld met een landspecifieke stekker voor aansluiting op het stroomnet van de faciliteit. Voor sommige landen, waaronder de Verenigde Staten, is de stekker stroomsterktespecifiek. Controleer vóór gebruik of de stroomspecificaties van uw faciliteit overeenkomen met die van de voedingskabel van de touchscreenconsole. De hardware van de mixer voor eenmalig gebruik (S.U.M.) van 2.000 I is niet geschikt voor circuits met een aardlekschakelaar (GFCI) omdat daarmee ongewenste uitschakeling kan optreden.

2.1.2 Voorbereiding van de hardware

De hardware wordt rechtstreeks door de fabrikant verzonden en wordt afgeleverd met verschillende veiligheidsmechanismen. Volg de onderstaande richtlijnen om de S.U.M. bij aankomst te configureren.



VOORZICHTIG: Alle procedures waarvoor de touchscreenconsole moet worden geopend, moeten worden uitgevoerd met een uitgeschakelde, vergrendelde hoofschakelaar en alle voedingsbronnen moeten van de touchscreenconsole zijn losgekoppeld. Voor de veiligheid van de gebruiker moet u de S.U.M.-hardware op zijn plek vastzetten door voorafgaand aan het onderhoud de zwenkwielen uit te schakelen.

2.2 Hardwaremontage

2.2.1 Hardware uit de kist halen

De S.U.M.-hardware wordt rechtstreeks door de fabrikant verzonden en geleverd in een krat. Volg de bijgeleverde uitpakinstructies en bewaar alle verpakkingsmateriaal voor eventueel toekomstig gebruik. Als u merkt dat er tijdens het vervoer schade is ontstaan, neemt u onmiddellijk contact op met uw salesmedewerker.

De S.U.M. wordt afgeleverd met de volgende items:

- Buitenste ondersteuningscontainer, met inbegrip van platform, tank, sondedeurplaten en touchscreenconsole (bevestigd).
- Powdertainer-arm-eenheid (optioneel)
- Aandrijfas

- Moment- en moersleutels
- Pompplanken (optioneel)
- Mand (optioneel)
- Kabelkapstok (optioneel)
- Kabelbeheerarm (optioneel)
- Communicatiekabels, met gereedschap om kabels aan te sluiten op de touchscreenconsole (afbeelding 2.1)
- Specificatiepakket van de apparatuur (ETP) op een USB-stick (apart verzonden)



Afbeelding 2.1. Gereedschap om kabels op de touchscreenconsole aan te sluiten.

Zie de Uitpakhandleiding voor HyPerforma S.U.M. met touchscreenconsole (DOC0061) voor meer informatie over het uitpakken van de S.U.M.

2.2.2 Motoreenheid voor S.U.M. van 2.000 I

Vanwege de omvang van de S.U.M.-eenheid van 2.000 I moet de mengmotor voor verzending worden verwijderd. Volg alleen voor S.U.M.-eenheden van 2.000 I de onderstaande stappen om de mengmotor op de buitenste ondersteuningscontainer te monteren. **Opmerking:** Voor de veiligheid dient de motor door twee personen op de tank te worden gemonteerd.

- 1. Haal het motor uit de doos.
- 2. Verwijder de bevestigde montageplaat van de veiligheidsdeksel van de bovenkant van de motor.
- 3. Zet de motor op het motorblokgedeelte van de ondersteuningsarm (afbeelding 2.2).



Motorblokgedeelte van ondersteuningsarm

Motorondersteuningsarm

Afbeelding 2.2. Motorondersteuningsarm voor S.U.M. van 2.000 I.

- 4. Terwijl de ene persoon de motor op zijn plek houdt, plaatst de tweede persoon de montageplaat van de veiligheidsdeksel op de bovenkant van de motor.
- 5. Sluit de veiligheidsdeksel van de motor en gebruik de pen (aan een kabel bevestigd) om deze vast te zetten.
- 6. Breng de vier bouten van 0,95 x 15,8 cm (3/8 x 6,25 inch) aan via de plaat, motor, en in het motorblokgedeelte van de ondersteuningsarm.
- 7. Draai alle vier de bouten aan.
- Voer de sensordraad van de veiligheidsdeksel naar de module met de wisselstroommotor achter de touchscreenconsole. Sluit de M12-stekker aan op het lege stopcontact op de onderkant van de module met de wisselstroommotor.
- 9. Voer de voedingskabel van de motor naar de module met de wisselstroommotor achter de touchscreenconsole. Sluit de rechthoekige stekker met zes pootjes aan op het lege stopcontact op de onderkant van de module van de wisselstroommotor en zet de verbinding vast met de hendelklem.

2.2.3 Voorbereiding van laadcellen

Wanneer S.U.M.-eenheden worden aangeschaft met in de fabriek geïnstalleerde laadcellen, worden de laadcellen in vergrendelde positie (omhoog gedraaid) verzonden om de apparatuur te beschermen (afbeelding 2.3).



Afbeelding 2.3. Details van laadcellen.

Na plaatsing van de apparatuur volgt u de onderstaande instructies om de laadcellen te ontgrendelen.

- 1. Verwijder de Delrin-sleepring als die aanwezig is.
- 2. Verwijder de drieklem.
- 3. Maak de vergrendelmoer los met het kleine uiteinde van het bijgeleverde gereedschap. Ga door totdat de moer strak tegen de basis of poot van de tank zit.
- 4. Herhaal dit proces voor elke laadcel totdat alle vergrendelmoeren los zijn van de vergrendelstangen.
- 5. Plaats de drieklem niet terug.

6. Na ontgrendeling moeten de laadcellen worden gekalibreerd in overeenstemming met uw bedrijfsbeleid. Zie paragraaf 4.5 voor instructies om de massa te kalibreren met de touchscreenconsole.

VOORZICHTIG: Verwijder de eenheid niet wanneer de laadcellen zijn ontgrendeld, en vooral niet wanneer de eenheid is gevuld. Hierdoor kunnen de laadcellen beschadigd raken.

Ontgrendelde laadcellen opnieuw vergrendelen:

- 1. Draai de vergrendelmoer met de hand vast op de stang.
- 2. Gebruik het meegeleverde gereedschap om de moer nog een kwartslag te draaien.

VOORZICHTIG: Om schade aan de laadcellen te voorkomen, mag u de moer niet te strak aandraaien.

- 3. Breng een standaard rvs drieklem van 38,1 mm (1,5 inch) aan rond de flenzen.
- 4. Herhaal dit proces voor alle laadcellen.

2.2.4 Extra items voor montage

 Als u sondes gaat gebruiken en optionele sondeclips (afbeelding 2.4) hebt besteld, bevestigt u de sondeclips aan de toegangsinkeping voor sondes op de klamp.



Afbeelding 2.4. Plastic sondeclip.

2. Plaats de aandrijfas, gereedschappen en lagerpot in de beugels van de S.U.M.-hardware-eenheid (afbeelding 2.5 en 2.6)



Afbeelding 2.5. Gereedschap in de beugel.



Afbeelding 2.6. Aandrijfas en lagerpot in de beugel.

3. Als uw systeem is voorzien van een watermantel, gebruikt u de aansluiting met drie klemmen om de in- en uitgangen van de watermantel aan te sluiten op onderkant van de buitenste ondersteuningscontainer (afbeelding 2.7 en 2.8).



Afbeelding 2.7. Aansluiting met drie klemmen op uitgang van watermantel verwijderen.



Afbeelding 2.8. Aansluiting op poort van watermantel bevestigen.

4. Als uw eenheid is voorzien van de optionele Powdertainer-arm, installeert u deze op de buitenste ondersteuningscontainer zoals hierna getoond (afbeelding 2.9).



Afbeelding 2.9. Powdertainer-arm installeren.

5. Als uw eenheid wordt geleverd met de (optionele) pompplanken (afbeelding 2.10), bevestigt u deze met een sleutel en vier bouten per plank aan de staaf. De (optionele) mand is vooraf aangebracht.



Afbeelding 2.10. Planken en mand bevestigd aan de S.U.M.

2.3 Hardwareconfiguratie

De S.U.M.-hardware mag alleen worden verplaatst over gladde oppervlakken. De S.U.M. moet leeg zijn en zijn losgekoppeld van alle stroom- en toevoerbronnen. Als uw systeem is voorzien van laadcellen, moeten deze worden vergrendeld voordat de S.U.M.eenheid wordt verplaatst. Volg de stappen hierna om de S.U.M.hardware voor gebruik te configureren.

 Verplaats de S.U.M. naar de gewenste plek en vergrendel de zwenkwielen om de eenheid te blokkeren. De S.U.M. van 2.000 l heeft geen wielen.

Opmerking: De eindgebruiker is ervoor verantwoordelijk alle bewegingen van het systeem te voorkomen en de geschiktheid van de structuur van het gebouw voor de betreffende apparatuur te beoordelen, om te verzekeren dat de verankering en de bevestiging van de apparatuur en het gebouw bestand zijn tegen seismische activiteit.

- 2. Controleer of de elektrische voedingen in de faciliteit voldoende zijn voor de stroomvereisten van de S.U.M. en bijbehorende componenten, zoals controllers of pompen.
- Als uw eenheid een watermantel heeft, sluit u de in- en uitgangen van de temperatuurcontrole-unit (TCU) aan op de watermantelpoorten van de S.U.M. (afbeelding 2.11 en 2.12). De ingang zit gewoonlijk links wanneer u naar de connectoren kijkt.



Afbeelding 2.11. De TCU aansluiten op de watermantel.



Afbeelding 2.12. Verbindingen tussen de S.U.M. en de TCU.

Opmerking: Raadpleeg de richtlijnen van de fabrikant van de TCU voor installatie- en bedieningsinstructies voor de TCU.

4. Als u een controller van derden gebruikt, zoekt u de te gebruiken digitale communicatiepoort op de achterkant van de touchscreenconsole op. Voordat u de aansluiting tot stand brengt, moet u controleren of de juiste T-stekker en terminator zijn aangebracht. Een kit met deze onderdelen is op bestelling verkrijgbaar. Raadpleeg paragraaf 8.5–Specificaties van accessoires en opties. U kunt een controller ook aansluiten met de verouderde poort op de achterkant van de touchscreenconsole.
3

Configuratie van de touchscreenconsole

Inhoudsopgave

- 3.1 De hardware van de touchscreenconsole configureren
- 3.2 Het startscherm van de touchscreenconsole configureren
- 3.3 Moduleoverzicht
- 3.4 Vereiste externe onderdelen voor modules
- 3.5 Modules configureren

3.1 De hardware van de touchscreenconsole configureren

Afbeelding 3.1 illustreert het achterpaneel van de touchscreenconsole. De verschillende poorten zijn voor uw gemak gelabeld.



Afbeelding 3.1. Poorten en aansluitingen op de achterkant van de touchscreenconsole.

- 1. Zet alle pompen die op de S.U.M. worden aangesloten op de bevestigde planken, indien aanwezig (afbeelding 3.2).
- 2. Sluit de communicatiekabels van de pomp(en) aan op de pomp-/ hulpuitgang op de achterkant van de touchscreenconsole en zet ze vast met het bijgeleverde gereedschap (afbeelding 3.3).



Afbeelding 3.2. Pompen op planken plaatsen.



- Als u het optionele kabelbeheersysteem gebruikt, leidt u de pompkabels door de clips die op de handgrepen op de achterkant van de S.U.M. zijn bevestigd. Sluit de pompen vervolgens aan op de juiste voedingsbron. **Opmerking:** Als u de pompen wilt kalibreren nadat ze zijn geconfigureerd, moet u eerst paragraaf 3.2–3.5 doornemen. Raadpleeg vervolgens paragraaf 4.1 – Pompen kalibreren.
- 4. Als u lucht wilt aansluiten op het S.U.M.-systeem, sluit dan een luchtleiding met een buitendiameter van 8 mm (5/16 inch) (afbeelding 3.4 — niet bijgeleverd) aan op de eerste persluchtpoort op de achterkant van de touchscreenconsole (afbeelding 3.1) en op de luchttoevoer van uw faciliteit. VOORZICHTIG: De luchtdruk naar de touchscreenconsole moet tussen 4,14–6,55 bar (60–95 psi) liggen.



Afbeelding 3.4. Luchtleiding aangesloten op de achterkant van de touchscreenconsole.

5. Controleer of de hoofdonderbreker en de hoofdschakelaar op 'uit' staan. De keuzeschakelaar moet wijzen naar de kleine cirkel links (afbeelding 3.5).



Afbeelding 3.5. Voedingsschakelaar in de 'uit'-stand.

 Verifieer dat de noodstop ('E-Stop') vrij is (uitgetrokken), door de knop naar rechts te draaien (afbeelding 3.6).
 Opmerking: De E-Stop verbreekt alle stroom naar het systeem. Een alarmzoemer gaat af wanneer de E-Stop wordt geactiveerd.



Afbeelding 3.6. De E-Stop rechtsom draaien om deze uit te schakelen.

7. Zorg dat de jumperstekker op de achterkant van de touchscreenconsole is aangesloten.

8. Sluit de voedingskabel van de motor aan op de motor (afbeelding 3.7).



Afbeelding 3.7. Aansluiting motorbesturingseenheid.

- Sluit alle elektrische stekkers aan op de stroomvoorziening van de vestiging. Raadpleeg de labels en de schema's van de hardware/ elektriciteit om te verzekeren dat de juiste elektrische spanning op de S.U.M. aangesloten is.
- 10. Gebruik het meegeleverde gereedschap (afbeelding 3.8) om alle communicatiekabels die u zult gebruiken (temperatuursensor, pH-sensor, druksensor, geleidbaarheidssensor, laadcellen enz.), aan te sluiten op de ingangspoorten op de achterzijde van de touchscreenconsole (afbeelding 3.9). **Opmerking:** Opdat de labels tijdens het maken van de aansluitingen zichtbaar blijven, raden we aan om met het aansluiten van de kabels op de touchscreenconsole onderaan te beginnen en naar boven toe te werken.



Afbeelding 3.8. Bijgeleverd gereedschap om communicatiekabels aan te sluiten.



Afbeelding 3.9. Communicatiekabels aansluiten op de touchscreenconsole.

11. Gebruik zo nodig de noodstroompoorten op de achterkant van de touchscreenconsole **Opmerking:** Noodstroom heeft dezelfde lijnspanning als de touchscreenconsole (120 of 240 V, afhankelijk van uw systeem). Verifieer voor het gebruik dat de hulponderbreker op de achterzijde van de touchscreenconsole op 'on' (aan) staat.

VOORZICHTIG: Alle vier de noodstroompoorten delen dezelfde onderbreker [8 ampère (A) stroom vanuit de touchscreenconsole]. Om doorslaan van de hulponderbreker te voorkomen mag u geen stroom van meer dan 8 A gebruiken voor de vier noodstroompoorten samen.

12. Wanneer u klaar bent om het systeem te gebruiken, schakelt u eerst de hoofdonderbreker in en daarna de hoofdschakelaar. Nadat de touchscreenconsole is opgestart, controleert u of de noodstop is uitgeschakeld en drukt u vervolgens op de blauwe knop 'stroom aan resetten' (afbeelding 3.10).



Afbeelding 3.10. Drukken op de blauwe resetknop om de stroom in te schakelen.

3.2 Het startscherm van de touchscreenconsole configureren

3.2.1 Kennismaking met de touchscreenconsole

Afbeelding 3.11 toont een voorbeeld van het startscherm van de touchscreenconsole. Zie tabel 3.1 voor meer informatie over de op het scherm genummerde items.



Afbeelding 3.11. Voorbeeld van het startscherm van de touchscreenconsole.

Tabel 3.1. Beschrijvingen van items op het startscherm van de touchscreenconsole.

Artikel	Naam	Beschrijving		
1	Beheerderspictogram	Als er een beheerdersprofiel is gemaakt en dit momenteel is aangemeld, dan is dit pictogram blauw. Als de beheerder niet is aangemeld, is het pictogram grijs.		
2	Alarmpictogram	Stuurt de gebruiker door naar het scherm Alarm Status (Alarmstatus), waar alarmen kunnen worden weergegeven/bevestigd.		
3	Instellingenpictogram	Stuurt de gebruiker door naar het scherm System Settings (Systeeminstellingen), waar de gebruiker de datum/tijd kan aanpassen, een beheerdersprofiel kan instellen, kan zien of er software-updates zijn enz.		
4	Uitklapbare menubalk	Toont de actieve modules. Door op de naam van de module in het menu te drukken, wordt de actieve module geopend.		
5	Geconfigureerde modules	Toont de modules die de gebruiker aan het startscherm heeft toegevoegd en geconfigureerd. Zie paragrafen 3.2.2 en 3.2.3 voor informatie over het toevoegen en configureren van modules.		
6	Grootte, datum en tijd van de S.U.M.	Toont de grootte van de momenteel gebruikte S.U.M. en de huidige datum en tijd.		
7	Paginanavigatie	Het startscherm bestaat uit drie pagina's, die onder aan het scherm zichtbaar zijn als drie stippen. De stip voor de actieve pagina is wit, de andere zijn gedimd. U kunt de volgende beschikbare pagina openen door op het startscherm van rechts naar links te vegen.		

3.2.2 Modules toevoegen aan het startscherm

De touchscreenconsole gebruikt 'modules' om de functies te identificeren die kunnen worden gebruikt. Paragraaf 3.2.3 bevat overzichtsinformatie over elke module. Volg de onderstaande stappen om de modules toe te voegen die u op het startscherm wilt gebruiken.

 Nieuwe gebruikers zien na het inschakelen van de touchscreenconsole een leeg scherm (afbeelding 3.12).
 Opmerking: Voordat u begint modules toe te voegen aan het startscherm, moet u zorgen dat de maateenheden, datum en tijd rechtsonder op het lege scherm juist zijn. Als de waarden onjuist zijn, raakt u het pictogram 'Instellingen' rechtsboven aan en selecteert u 'Instrument Settings' (Instrumentinstellingen) om de waarden in te stellen.



Afbeelding 3.12. Bij het opstarten verschijnt een leeg scherm.

- 2. Wanneer de maateenheden, datum en tijd juist op het scherm worden weergegeven, raakt u 'Home screen' (Startscherm) aan.
- In het geopende scherm Select Auxiliary Outputs or Pumps (Hulpuitgangen of pompen selecteren) kan de gebruiker een combinatie van vier pompen en/of hulpuitgangen selecteren (afbeelding 3.13). Raak de pompen en/of hulpuitgangen aan die u wilt toevoegen aan uw startscherm. Druk opnieuw op de knop om de selectie van een pomp of hulpuitgang ongedaan te maken. Wanneer u gereed bent, raakt u 'Next' (Volgende) aan.



Afbeelding 3.13. Pompen en/of hulpuitgangen selecteren.

4. In het geopende scherm Select Modules (Modules selecteren) kan de gebruiker andere modules selecteren (afbeelding 3.14). Raak de modules aan die u wilt toevoegen aan het startscherm. Druk opnieuw op de knop om de selectie van een module ongedaan te maken. Wanneer u gereed bent, raakt u 'Next' (Volgende) aan.



Afbeelding 3.14. Modules selecteren.

 In het geopende scherm Home Screen Confirmation (Startscherm bevestigen) worden de pompen/hulpuitgangen en modules weergegeven die u hebt geselecteerd (afbeelding 3.15). Klik op 'Confirm' (Bevestigen) om ze aan uw startscherm toe te voegen. Het volgende geopende scherm ziet er vergelijkbaar uit, maar bevat de modules die aan het startscherm zijn toegevoegd. Raak 'Open Home Screen' (Startscherm openen) aan.

€	Workspace Confirmation		
	Confirm that all the modules for	or this workspace are correct	
	Agitation AuxIn2 BPC pressure Fill Pump1 Pump4 Timer	AuxIn1 AuxOut3 Conductivity Mass Pump2 Temperature pH	
		Cancel Confirm	

Afbeelding 3.15. Bevestigingsscherm van startscherm.

 Het geconfigureerde startscherm wordt geopend met de door u geselecteerde modules. Al naar gelang het aantal toegevoegde modules kan uw startscherm uit meerdere pagina's bestaan. Veeg van rechts naar links om andere pagina's weer te geven. Zie de volgende paragraaf, 3.3 – Moduleoverzicht, voor algemene informatie over de functionaliteit van elke module en instructies om modules te configureren.

3.2.3 Een beheerdersprofiel maken en gebruiken

Met een beheerdersprofiel kan een beheerder de toegang van andere gebruikers tot bepaalde instellingen en functies van de touchscreenconsole beperken. Tabel 3.2 toont de beschikbare functies voor 'alle gebruikers' en 'alleen beheerders'.

Instelling of functie	Alle gebruikers	Alleen beheerders	
pH-instellingen	Calibration (Kalibratie), Temperature compensation (Temperatuurcompensatie), Select active sensor (Actieve sensor selecteren), Reset calibration (Kalibratie resetten), Module configuration (Moduleconfiguratie) en Sensors available (Beschikbare sensoren)	Alarmen High (Hoog) en Low (Laag), Alarmen High High (Hoog-hoog) en Low Low (Laag-laag), Alarm designation (Alarmaanduiding) (pompen) en Diagnostics (Diagnose)	
Systeeminstellingen	Export sensor logs (Sensorlogboeken exporteren)	About instrument (Info instrument), Maintenance and services (Onderhoud en services), Instrument settings (Instrumentinstellingen), Alarms and interlocks (Alarmen en interlocks) en End User License Agreement (Gebruiksrechtovereenkomst)	
BPC-drukinstellingen	Units (Eenheden), Calibration (Kalibratie)	Alarmen High (Hoog) en Low (Laag), Alarmen High High (Hoog-hoog) en Low Low (Laag-laag), Alarm designation (Alarmaanduiding) (pompen) en Module Configuration (Module configureren)	
Roer-instellingen	N.v.t.	Alarmen High (Hoog) en Low (Laag), Alarmen High High (Hoog-hoog) en Low Low (Laag-laag), Set point limits (Limieten instelpunten), Calibration (Kalibratie) en Module configuration (Module configureren)	
Temperatuurinstellingen	Select active sensor (Actieve sensor selecteren), Calibration (Kalibratie)	Alarmen High (Hoog) en Low (Laag), Alarmen High High (Hoog-hoog) en Low Low (Laag-laag), TCU set point limits (Instelpuntlimieten TCU) en Module configuration (Module configureren)	
Timerinstellingen	Notes (Opmerkingen), Reset (Resetten)	N.v.t.	
Massa-instellingen	Fluid density (Vloeistofdichtheid), Tare/zero (Tarreren/nulstelling)	Alarmen High (Hoog) en Low (Laag), Alarmen High High (Hoog-hoog) en Low Low (Laag-laag), Alarm designation (Alarmaanduiding) (pompen), Alarm designation (Alarmaanduiding) (roeren), Alarm designation (Alarmaanduiding) (knijpventielen), Calibration (Kalibratie) en Module configuration (Module configureren)	
Pompinstellingen	Units and fluid density (Eenheden en vloeistofdichtheid), Pump calibration (Pompkalibratie), Add bolus (Bolus toevoegen) en Prime pump (Pomp primen)	Diagnostics (Diagnose), RPM scaling (RPM-schaling), RPM limits (RPM-limieten) en Module configuration (Module configureren)	
Vloeistofdrukinstellingen	Units (Eenheden), Display options (Weergaveopties) en Calibration (Kalibratie)	Low and High alarms (Alarmen Laag en Hoog), Low Low and High High Alarms (Alarmen (Low Low) Laag-laag) en High High (Hoog-hoog)), Alarm designation (Alarmaanduiding) (pompen), Alarm designation (Alarmaanduiding) (roeren), Alarm designation (Alarmaanduiding) (knijpventielen) en Module configuration (Module configureren)	

Tabel 3.2. Toegankelijkheid van instellingen/functies voor alle gebruikers en alleen behee	rders.
--	--------

Om een beheerdersprofiel in te stellen voert u de volgende stappen uit.

- 1. Open 'System Settings' (Systeeminstellingen) door het pictogram 'Settings' (Instellingen) rechtsboven in het startscherm aan te raken. Raak 'Administrator' (Beheerder) aan.
- Voer in het geopende scherm de pincode in die u als beheerder wilt gebruiken. Voer de pincode opnieuw in bij 'Confirm PIN' (Pincode bevestigen). **Opmerking:** Als u de door u ingevoerde pincode wilt weergeven, raakt u het vak 'Show PIN' (Pincode weergeven) aan.
- Wanneer u klaar bent, raakt u 'Create profile' (Profiel maken) aan. Er verschijnt een bevestiging dat uw beheerdersprofiel is gemaakt. Raak 'Done' (Gereed) aan.
- Nadat u een beheerdersprofiel hebt gemaakt, kunt u zich met de optie 'Administrator' (Beheerder) onder 'System settings' (Systeeminstellingen) bij het profiel af- en aanmelden (afbeelding 3.16).



Afbeelding 3.16. Scherm Administrator profile (Beheerdersprofiel).

Selecteer 'Actions' (Acties) voor meer beheerdersopties. De volgende opties zijn beschikbaar:

- Reset PIN (PIN terugzetten): de PIN van het beheerdersprofiel wijzigen.
- Enable auto sign out (Automatisch afmelden inschakelen): toestaan dat de touchscreenconsole de beheerder automatisch afmeldt na een door de gebruikers opgegeven tijdsduur.

• Control admin PIN (Beheerders-PIN controleren): het beheerdersprofiel activeren of deactiveren. **Opmerking:** Als u 'Disable' selecteert, wordt het beheerdersprofiel uit de touchscreenconsole verwijderd. Als u het profiel wilt behouden maar nu niet wilt gebruiken, meldt u zich in plaats daarvan af bij het profiel.

3.3 Moduleoverzicht

Met elk van de hierna vermelde touchscreenconsole-modules kunnen gebruikers verschillende functies bewaken en instellen. Sommige modules vereisen, of kunnen optioneel worden gebruikt met, hulpcomponenten zoals sensoren of pompen. Veel modules zijn ook 'optioneel afhankelijk' van hulpcomponenten of andere modules. Dit betekent dat gebruikers desgewenst binnen modules interlocks kunnen instellen die hun functie afhankelijk van bepaalde voorwaarden wijzigen, zoals het uitschakelen van een pomp wanneer de BPCluchtdruk een bepaalde PSI-waarde bereikt. Zie paragraaf 3.5.1 voor meer informatie om interlocks in te stellen.

- Agitation (Roeren) Deze module wordt gebruikt om de mengsnelheid van de S.U.M. te bewaken en regelen. Deze snelheid wordt weergegeven in omwentelingen per minuut (RPM). De module Agitation (Roeren) is optioneel afhankelijk van de temperatuurcontrole-unit (TCU), als er een TCU wordt gebruikt. Dit betekent dat de touchscreenconsole de TCU kan pauzeren of stoppen als de roersnelheid boven of onder de gespecificeerde instelpunten komt.
- Mass (Massa) deze module heeft laadcellen nodig om het totale gewicht van het S.U.M-systeem te lezen. De module Mass (Massa) is optioneel afhankelijk van het volgende:
 - Knijpventiel voor vullen: Het knijpventiel sluit als de massaaflezing een gespecificeerd instelpunt overschrijdt, om overlopen te voorkomen.
 - Knijpventiel voor aftappen: Het knijpventiel sluit als de massaaflezing onder een gespecificeerd instelpunt komt, om teveel aftappen te voorkomen.
 - Pompen: De pompen worden uitgeschakeld als de massaaflezing boven of onder een gespecificeerd instelpunt komt omdat de knijpventielen voor vullen of aftappen sluiten.
 Dit voorkomt overdruk in BPC-lijnen en barsten daarvan.
 - Module Agitation (Roeren): Het roeren stopt als de massaaflezing onder een gespecificeerd instelpunt komt, om potentiële schade voor mengen in een lege BPC te voorkomen.

- 3. BPC Pressure (BPC-druk) Deze module kan alleen worden gebruikt met closed-top BPC's die zijn voorzien van druksensoren, en in faciliteiten waarin een luchtbron beschikbaar is. BPC Pressure (BPC-druk) wordt gebruikt om de BPC te vullen met een gespecificeerde hoeveelheid lucht en om de druk in de hoofdruimte van de BPC te bewaken. De functie is optioneel afhankelijk van het knijpventiel voor vullen en de pompen: het knijpventiel sluit en de pomp stopt als de psi boven een gespecificeerd instelpunt komt.
- 4. Pumps (Pompen) Deze module kan alleen worden gebruikt als een of meer pompen ingesteld zijn. De module Pumps (Pompen) kan worden gebruikt om de BPC te vullen met vloeistof, zuur-, base- of andere bufferoplossingen te leveren en om vloeistof af te tappen. Het gebruik van knijpventielen is optioneel, maar bij gebrek daaraan moeten gebruikers BPC-lijnen handmatig openen en sluiten met knijpventielen.
- 5. Fill (Vullen) Deze module kan alleen worden gebruikt als uw systeem is uitgerust met laadcellen en er een of meer pompen zijn geconfigureerd. De module Fill (Vullen) wordt gebruikt om de BPC voor het mengen met vloeistof te vullen. Het gebruik van knijpventielen is optioneel, maar zonder ventielen moeten gebruikers de BPC vullen door handmatig een knijpventiel op een BPC-leiding te openen.
- 6. Harvest (Aftappen) Deze module kan alleen worden gebruikt als uw systeem is uitgerust met laadcellen en er een pomp is geconfigureerd. De module Harvest (Aftappen) wordt gebruikt om vloeistof uit de BPC af te tappen na voltooiing van het mengproces. Het gebruik van knijpventielen is optioneel, maar zonder ventielen moeten gebruikers de BPC aftappen door handmatig een knijpventiel op een BPC-leiding te openen.
- 7. pH Deze module kan niet worden gebruikt zonder een of twee pH-sensoren. De pH-module kan worden gebruikt om de pHniveaus tijdens het mengproces te bewaken en aan te passen. De mogelijkheid om pH-niveaus aan te passen is echter afhankelijk van het gebruik van pompen (bijv. één om een zuuroplossing af te leveren en een andere voor een base-oplossing). De module is optioneel afhankelijk van pompen omdat de pomp kan worden gepauzeerd of gestopt als de pH-waarde boven of onder gespecificeerde instelpunten komt.

- 8. Conductivity (Geleidbaarheid) Deze module kan niet worden gebruikt zonder een of twee geleidbaarheidssensoren. De module Conductivity (Geleidbaarheid) kan worden gebruikt om de geleidbaarheidsniveaus tijdens het mengproces te bewaken en aan te passen. De mogelijkheid om geleidbaarheidsniveaus aan te passen is echter afhankelijk van het gebruik van een pomp om een bufferoplossing af te leveren. De module is optioneel afhankelijk van pompen omdat de pomp kan worden gepauzeerd of gestopt als de geleidbaarheidswaarde boven of onder gespecificeerde instelpunten komt.
- 9. Temperature (Temperatuur) Deze module kan niet worden gebruikt zonder een of twee temperatuursensoren. De module Temperature (Temperatuur) kan worden gebruikt om de temperatuur te bewaken en aan te passen. De mogelijkheid om de temperatuur aan te passen is echter afhankelijk van het gebruik van een TCU. De module Temperature (Temperatuur) is optioneel afhankelijk van de TCU omdat de TCU stand-by kan worden gezet als de temperatuur boven of onder gespecificeerde instelpunten komt.
- Timer De module Timer bestaat uit een timer die bij nul begint en oploopt om te volgen hoe lang een oplossing wordt gemengd. De module heeft start- en stopfuncties.
- 11. Liquid Pressure (Vloeistofdruk) deze module kan niet worden gebruikt zonder een of twee vloeistofdruksensoren in de slangenset van de BPC (voor en/of na het filter). De monitor Liquid Pressure (Vloeistofdruk) wordt gebruikt om het vloeistofdrukniveau tijdens het aftappen te bewaken om te zorgen dat het filter niet verstopt raakt. De module Liquid Pressure (Vloeistofdruk) is optioneel afhankelijk van beide pompen en de knijpventielen voor aftappen omdat het knijpventiel sluit en de pomp stopt als de vloeistof boven een gespecificeerd instelpunt komt.
- 12. Auxiliary Output (Hulpuitgang) deze module kan worden gebruikt om indien nodig een andere bestaande module, zoals Temperature (Temperatuur) te dupliceren. De module Auxiliary Output (Hulpuitgang) gebruikt dezelfde verbindingspoort (op de achterzijde van de touchscreenconsole) als de pompen, zodat gebruikers een maximale combinatie van vier pompen en/of hulpuitgangen kunt hebben.
- Auxiliary Inputs (Hulpingangen) deze module kan worden gebruikt om sensoren en zenders te monitoren die een 4-20 mAsignaal gebruiken, zoals externe weegschalen.

3.4 Vereiste externe onderdelen voor modules

Het hierna afgebeelde stroomdiagram (afbeelding 3.17) illustreert de vereiste externe onderdelen voor verschillende modules.



Afbeelding 3.17. Stroomdiagram met verplichte externe componenten voor elke module.

3.5 Modules configureren

Elke moduleconfiguratie vereist dat de gebruiker bepaalde velden (zoals instelpunten) invult en bevat optionele velden (zoals alarmen) voor aanpassing. De volgende paragrafen bevatten informatie om interlocks, alarmen en individuele modules in te stellen.

3.5.1 Interlocks en alarmen instellen

Gebruikers kunnen de functionaliteit van hun touchscreenconsole aanpassen door interlocks en alarmen op individuele modules in te stellen.

- Door het instellen van 'interlocks' binnen een module ontstaan afhankelijkheden van andere modules of hulpcomponenten. Gebruikers kunnen bijvoorbeeld interlocks instellen in de module Mass (Massa) waarmee roeren wordt uitgeschakeld als de massa van de S.U.M. een bepaald gewicht bereikt. Deze interlock voorkomt potentiële schade die kan ontstaan door het mengen in een lege BPC of het mengen met een snelheid buiten de aanbevolen specificaties voor veilig gebruik van de aandrijfas.
- Alarmen zijn ook optioneel op modules van de touchscreenconsole. Gebruikers kunnen 'High' (Hoog) en 'Low' (Laag) alarmen instellen, of 'High High' (Hoog-hoog) en 'Low Low' (Laag-laag) alarmen.
 - De alarmen High (Hoog) en Low (Laag) waarschuwen de gebruiker alleen visueel dat de gespecificeerde instelpunten zijn bereikt. Wanneer de alarmen High (Hoog) of Low (Laag) worden geactiveerd, worden het alarmpictogram en de betreffende module op het startscherm geel weergegeven, maar worden de processen voortgezet.
 - De alarmen High High (Hoog-hoog) en Low Low (Laaglaag) waarschuwen de gebruiker visueel en hoorbaar dat de gespecificeerde instelpunten zijn bereikt. Wanneer de alarmen High High (Hoog-hoog) of Low Low (Laag-laag) worden geactiveerd, worden het alarmpictogram en de betreffende module rood weergegeven op het startscherm, klinkt er een zoemer en worden de processen gepauzeerd of gestopt (afhankelijk van de instellingen van de gebruiker).

Interlock- en alarminstellingen voor individuele modules zijn toegankelijk door de module op het startscherm aan te raken. Het scherm 'Configuration' (Configuratie) van elke module bevat velden om beide opties in te stellen. Zie voor meer informatie de instructies om modules te configureren in de volgende paragrafen.

3.5.2 De module Agitation (Roeren) configureren

Gebruik de volgende informatie en stappen om de module Agitation (Roeren) te configureren. Elk van de onderstaande items wordt geïllustreerd in afbeelding 3.18. **Opmerking:** U kunt 'Done' (Gereed) onder aan het scherm Configuration (Configuratie) te allen tijde aanraken tijdens de configuratie van de module. Uw invoer wordt opgeslagen en u kunt doorgaan waar u was gebleven wanneer u de moduleconfiguratie opnieuw opent.

Verplichte velden

• Set point limits (Limieten instelpunten): Voer de maximum- en minimumsnelheden (in RPM) in die u tijdens het mengproces wilt gebruiken.

Optionele velden

- Low and High alarms (Alarmen Laag en Hoog): Voer de RPM-waarden in waarmee de alarmen High (Hoog) en Low (Laag) worden geactiveerd. Deze alarmen hebben geen invloed op lopende processen, maar waarschuwen gebruikers dat bepaalde omstandigheden buiten de gespecificeerde gebruiksparameters vallen.
- Low Low and High High alarms (Alarmen Laag-laag en Hooghoog): Vul de velden voor de interlockvergelijking en hysterese in om de voorwaarden in te stellen waarmee de alarmen High High (Hoog-hoog) en Low Low (Laag-laag) worden geactiveerd. Bij activering van deze alarmen klinkt een zoemer en worden lopende processen gepauzeerd of gestopt, afhankelijk van de gebruikersinstellingen. De interlockvergelijking geeft aan onder welke omstandigheden het alarm wordt geactiveerd, terwijl het hysteresegedeelte aangeeft onder welke omstandigheden processen worden hervat. Raak elk van de getoonde velden aan om de vereiste waarden voor de alarmen High High (Hoog-hoog) en/of Low Low (Laag-laag) in te voeren of te selecteren.
- Calibration (Kalibratie) (alleen voor systemen met wisselstroommotoren): Stel éénpunts roerkalibratie in om de nauwkeurigheid van de roeraflezingen te bepalen. Zie hoofdstuk 4 voor meer informatie over kalibratieprocedures.
- Zoek de module Agitation (Roeren) op het startscherm van de touchscreenconsole op en raak 'Configure' (Configureren) aan. Het scherm Agitation Configuration (Roeren configureren) verschijnt (afbeelding 3.18). Raak het veld 'Set point limits' (Limieten instelpunten) aan.

Agitation Module Configuration					
Set Point Limits:	20RPM - 356RPM	•			
Low and High Alarms:	Lo Disabled-Hi Disab	~			
Low Low and High High Alarms:	LoLo Disabled-HiHi	~			
Calibration:	Calibrate	~			
		Done			

Afbeelding 3.18. Scherm Agitation Configuration (Roeren configureren). Gebruikershandleiding HyPerforma S.U.M. met touchscreenconsole | 49

- Voer in het geopende scherm de gewenste boven- en ondergrenzen voor de roer-instelpunten (in RPM) in. Raak 'Next' (Volgende) aan.
- Het scherm Agitation Configuration (Roeren configureren) wordt opnieuw geopend. De 'Max' en 'Min' waarden die u hebt opgegeven, moeten zichtbaar zijn in het veld naast 'Set point limits' (Limieten instelpunten). Als u de roer-alarmen High (Hoog)/ Low (Laag) wilt instellen voor roeren in relatie tot temperatuur, raakt u het veld 'Low and High alarms' (Alarmen Laag en Hoog) aan.
- 4. Het geopende scherm bevat aparte gedeeltes voor de alarmen High (Hoog) en Low (Laag) (afbeelding 3.19). Als u het alarm Low (Laag) wilt configureren, schuift u de knop On/Off (Aan/Off) naar de kant 'On' (Aan). Voer in het veld bij 'High Alarm' (Hoog alarm) de temperatuur [in graden Celsius (°C)] in waarbij het alarm High (Hoog) wordt geactiveerd.



Afbeelding 3.19. Scherm High and Low alarms (Alarmen hoog en laag).

- Als u het alarm Low (Laag) wilt configureren, schuift u de knop On/Off (Aan/Off) naar de kant 'On' (Aan). Voer in het veld bij 'Low Alarm' (Laag alarm) de temperatuur [in graden Celsius (°C)] in waarbij het alarm Low (Laag) wordt geactiveerd.
- 6. Nadat u de waarden voor de alarmen High (Hoog) en Low (Laag) hebt ingevoerd, raakt u 'Done' (Gereed) aan.
- 7. In het scherm Agitation Configuration (Roeren configureren) worden nu de ingevoerde waarden voor de alarmen High (Hoog) en Low (Laag) weergegeven. Als u de alarmen High High (Hoog-

hoog)/Low Low (Laag-laag) wilt instellen voor roeren in relatie tot temperatuur, raakt u het veld 'Low Low and High High alarms' (Alarmen Laag-laag en Hoog-hoog) aan.

 Het geopende scherm (afbeelding 3.20) bevat aparte gedeeltes voor de alarmen High High (Hoog-hoog) en Low Low (Laag-laag). Schuif de knoppen On/Off (Aan/uit) in elk gedeelte naar de kant 'On' (Aan) om de alarmen in te schakelen.

٠	Agitati	Agitation High High and Low Low Alarms			
	High High Alarm	Off			
	Interlock:	Agitation>0RPM then Stop TCU			
	Hysteresis:	Resume if Agitation<0RPM for 0s			
	Low Low Alarm	Off O			
	Interlock:	Agitation<0RPM then Stop TCU			
	Hysteresis:	Resume if Agitation>0RPM for 0s			
		Cancel Done			

Afbeelding 3.20. Scherm High High and Low Low Alarms (Alarmen Hoog-hoog en Laag-laag).

- Voor de alarmen High High (Hoog-hoog) en Low Low (Laaglaag) worden er velden met het label 'Interlock' en 'Hysteresis' (Hysterese) weergegeven. Raak 'Interlock' in een van de alarmvelden aan om de initiële voorwaarde in te stellen waarbij het alarm wordt geactiveerd, en aan te geven welke actie na activering moet worden uitgevoerd.
- 10. In het geopende scherm worden de interlock- en hysteresegedeeltes getoond. In het gedeelte 'TCU Interlock Equation' (TCU-interlockvergelijking) voert u waarden in bij Control Set Point (Controle-instelpunt) (in RPM) en selecteert u een waarde bij Action Taken (Ondernomen actie) (Pause (Pauzeren) of Stop (Stoppen)).
- 11. Het gedeelte 'Hysteresis' (Hysterese) (onder het gedeelte Interlock) bepaalt hoeveel RPM de S.U.M. moet bereiken (en hoe lang die RPM moet worden aangehouden, in seconden) voordat de TCU wordt hervat. Voer een waarde in bij Set Point (Instelpunt) (in RPM) en Seconds (Seconden). Wanneer u gereed bent, raakt u 'Done' (Gereed) aan.

- Wanneer u de interlock en hysterese voor het alarm High High (Hoog-hoog) en Low Low (Laag-laag) hebt ingesteld, raakt u 'Done' (Gereed) aan in het scherm 'Agitation High High and Low Low Alarms' (Alarmen Hoog-hoog en Laag-laag voor roeren).
- 13. Als u het systeem gebruikt met een wisselstroommotor en een roerkalibratie wilt instellen, raakt u het veld 'Calibration' (Kalibratie) aan. Volg de kalibratie-instructies uit hoofdstuk 4 Kalibratieprocedures. U kunt de kalibratie ook overslaan en 'Done' (Gereed) aanraken om terug te gaan naar het startscherm.

3.5.3 De module Mass (Massa) configureren

Gebruik de volgende informatie en stappen om de module Mass (Massa) te configureren. Elk van de onderstaande items wordt geïllustreerd in afbeelding 3.21. **Opmerking:** U kunt 'Done' (Gereed) onder aan het scherm Configuration (Configuratie) te allen tijde aanraken tijdens de configuratie van de module. Uw invoer wordt opgeslagen en u kunt doorgaan waar u was gebleven wanneer u de moduleconfiguratie opnieuw opent.

Verplichte velden

• Units (Eenheden): Voer in welke eenheden u wilt gebruiken wanneer u de S.U.M. weegt [gram (g) of kilogram (kg)].

Optionele velden

- Low and High alarms (Alarmen Laag en Hoog): Voer de massawaarden (g of kg) in waarmee de alarmen High (Hoog) en Low (Laag) worden geactiveerd. Deze alarmen hebben geen invloed op lopende processen, maar waarschuwen gebruikers dat bepaalde omstandigheden buiten de gespecificeerde gebruiksparameters vallen.
- Low Low and High High alarms (Alarmen Laag-laag en Hooghoog): Vul de velden voor de interlockvergelijking en hysterese in om de voorwaarden in te stellen waarmee de alarmen High High (Hoog-hoog) en Low Low (Laag-laag) worden geactiveerd. Bij activering van deze alarmen klinkt een zoemer en worden lopende processen gepauzeerd of gestopt, afhankelijk van de gebruikersinstellingen. De interlockvergelijking geeft aan onder welke omstandigheden het alarm wordt geactiveerd, terwijl het hysteresegedeelte aangeeft onder welke omstandigheden processen worden hervat. Raak elk van de getoonde velden aan om de vereiste waarden voor de alarmen High High (Hoog-hoog) en/of Low Low (Laag-laag) in te voeren of te selecteren.

- Alarm Designation (pumps, pinch valves, and agitation) (Alarmaanduiding; pompen, knijpventielen en roeren): Nadat u de massa-alarmen High High (Hoog-hoog) en/of Low Low (Laaglaag) hebt ingesteld, gebruikt u deze drie velden om te selecteren op welke pompen u deze alarmen wilt toepassen. Bijvoorbeeld voor Pomp1 schakelt u het alarm Low Low (Laag-laag) in en het alarm High High (Hoog-hoog) uit.
- Calibration (Kalibratie): Stel éénpunts massakalibratie in om de nauwkeurigheid van de massalezingen van laadcellen te bepalen. Zie hoofdstuk 4 voor meer informatie over kalibratieprocedures.
- Zoek de module 'Mass' (Massa) op het startscherm van de touchscreenconsole op en raak vervolgens 'Configure' (Configureren) aan. Het scherm Mass Configuration (Massa configureren) verschijnt (afbeelding 3.21). Raak het veld 'Units' (Eenheden) aan.



Afbeelding 3.21. Scherm Mass Configuration (Massa configureren).

- 2. Selecteer gram of kilogram in het geopende scherm. Raak 'Next' (Volgende) aan.
- Het scherm Mass Configuration (Massa configureren) wordt opnieuw geopend. De geselecteerde eenheden worden weergegeven in het veld naast 'Units' (Eenheden). Als u de alarmen High (Hoog)/Low (Laag) wilt instellen voor massa, raakt u het veld 'Low and High alarms' (Alarmen Laag en Hoog) aan.
- 4. Het geopende scherm bevat aparte gedeeltes voor de alarmen High (Hoog) en Low (Laag) (afbeelding 3.22). Als u het alarm High (Hoog) wilt configureren, schuift u de knop On/Off (Aan/Uit) naar de kant 'On' (Aan). Voer in het veld bij 'High Alarm' (Hoog alarm) de massawaarde in waarbij het alarm High (Hoog) wordt geactiveerd.

۲		Mass High and I	_ow Alarms		
Would you like to set the High and Low alarms?					
			High Alarm		
	Off	On	0.0		
			kg		
			Low Alarm		
	Off	On	0.0		
			kg		
			Cancel	Done	

Afbeelding 3.22. Scherm High and Low alarms (Alarmen hoog en laag).

- Als u het alarm Low (Laag) wilt configureren, schuift u de knop On/Off (Aan/Off) naar de kant 'On' (Aan). Voer in het veld bij 'Low Alarm' (Laag alarm) de massawaarde in waarbij het alarm Low (Laag) wordt geactiveerd.
- 6. Nadat u de waarden voor de alarmen High (Hoog) en Low (Laag) hebt ingevoerd, raakt u 'Done' (Gereed) aan.
- 7. In het scherm Mass Configuration (Massa configureren) worden nu de ingevoerde waarden voor de alarmen High (Hoog) en Low (Laag) weergegeven. Als u de alarmen High High (Hoog-hoog)/Low Low (Laag-laag) wilt instellen voor massa (in verband met pompen, knijpventielen of roeren), raakt u het veld 'Low Low and High High alarms' (Alarmen Laag-laag en Hoog-hoog) aan.
- Het geopende scherm (afbeelding 3.23) bevat aparte gedeeltes voor de alarmen High High (Hoog-hoog) en Low Low (Laag-laag). Schuif de knoppen On/Off (Aan/uit) in elk gedeelte naar de kant 'On' (Aan) om de alarmen in te schakelen.

۲	Mass1 High High and Low Low Alarms		
	High High Alarm	Off Off	
	Interlock:	Mass1>0.0kg then Stop	
	Hysteresis:	Resume if Mass1<0.0kg for 0s	
	Low Low Alarm	Off Off	
	Interlock:	Mass1<0.0kg then Stop	
	Hysteresis:	Resume if Mass1>0.0kg for 0s	
		Cancel Done	

Afbeelding 3.23. Scherm High High and Low Low Alarms (Alarmen Hoog-hoog en Laag-laag).

- 9. Voor de alarmen High High (Hoog-hoog) en Low Low (Laaglaag) worden er velden met het label 'Interlock' en 'Hysteresis' (Hysterese) weergegeven. Raak 'Interlock' in een van de alarmvelden aan om de initiële voorwaarde in te stellen waarbij het alarm wordt geactiveerd, of 'Hysteresis' (Hysterese) om aan te geven welke actie na activering moet worden uitgevoerd.
- 10. In het geopende scherm worden de interlock- en hysteresegedeeltes getoond. In het gedeelte 'Agitation Interlock Equation' (Interlockvergelijking voor roeren) voert waarden in bij Control Set Point (Controle-instelpunt) (in g of kg) en selecteert u een waarde bij Action Taken (Ondernomen actie) (Pause (Pauzeren) of Stop (Stoppen)).
- 11. Het gedeelte 'Hysteresis' (Hysterese) (onder het gedeelte Interlock) bepaalt de massa die de S.U.M. moet bereiken (en hoe lang die massa moet worden aangehouden, in seconden) voordat het roeren, de pompen of de knijpventielen opnieuw worden geactiveerd. Voer een waarde in bij Set Point (Instelpunt) (in g of kg) en Seconds (Seconden). Wanneer u gereed bent, raakt u 'Done' (Gereed) aan.
- Nadat u de interlock- en hystereseparameters voor de alarmen Low Low (Laag-laag) en High High (Hoog-hoog) hebt ingesteld, raakt u 'Done' (Gereed) aan. Het scherm Mass Configuration (Massa configureren) wordt opnieuw geopend.
- 13. Als u ingestelde alarmen voor roeren of specifieke pompen/ knijpventielen wilt aanduiden, raakt u het bijbehorende veld 'Alarm Designation' (Alarmaanduiding) in het scherm Mass Configuration (Massaconfiguratie) aan. **Opmerking:** Wanneer een pomp of

knijpventiel is ingeschakeld, kan de massa worden aangepast nadat het alarm is geactiveerd. Bij inschakeling van een alarm voor roeren wordt het roeren in- of uitgeschakeld nadat het alarm is geactiveerd.

14. In het geopende scherm (afbeelding 3.24) ziet u de interlock- en hystereseparameters die u hebt opgegeven voor de alarmen High High (Hoog-hoog) en Low Low (Laag-laag). Eronder worden het roeren, de pompen of de knijpventielen weergegeven die zijn ingesteld met het systeem, afhankelijk van het veld Alarm Designation (Alarmaanduiding) dat u in het scherm Mass Configuration (Massaconfiguratie) hebt geselecteerd. Raak een knop voor een pomp, knijpventiel of roeren aan om de alarmen High High (Hoog-hoog) en Low Low (Laag-laag) in of uit te schakelen. **Opmerking:** Afbeelding 3.24 toont de opties voor de selectie 'Alarm Designation (pumps)' (Alarmaanduiding (pompen)).



Afbeelding 3.24. Scherm Mass Alarm Designation (Pumps) Alarmaanduiding Massa (pompen).

- 15. Schuif op het geopende scherm de knop Enable/Disable (In-/ uitschakelen) naar de kant Enable (Inschakelen) voor de alarmen High High (Hoog-hoog) en/of Low Low (Laag-laag) om het of de alarmen in te schakelen, voor roeren of voor de geselecteerde pomp of het geselecteerde knijpventiel. Wanneer u gereed bent, raakt u de knop 'Done' (Gereed) aan.
- 16. Als u een massakalibratie wilt instellen, raakt u het veld 'Calibration' (Kalibratie) aan. Volg de kalibratie-instructies uit hoofdstuk 4 — Kalibratieprocedures. U kunt de kalibratie ook overslaan en 'Done' (Gereed) aanraken om terug te gaan naar het startscherm.

3.5.4 De module Timer configureren

Omdat de module Timer alleen de functies 'Start' (Starten) en 'Stop' (Stoppen) bevat, hoeft deze module niet te worden geconfigureerd.

3.5.5 De module BPC Pressure (BPC-druk) configureren

Gebruik de volgende informatie en stappen om de module BPC Pressure (BPC-druk) te configureren. Elk van de onderstaande velden wordt geïllustreerd in afbeelding 3.25. **Opmerking:** U kunt 'Done' (Gereed) onder aan het scherm Configuration (Configuratie) te allen tijde aanraken tijdens de configuratie van de module. Uw invoer wordt opgeslagen en u kunt doorgaan waar u was gebleven wanneer u de moduleconfiguratie opnieuw opent.

Verplichte velden

• Units (Eenheden): Geef de eenheden op die u wilt gebruiken wanneer u de BPC vult met lucht [millibar (mbar) of psi].

Optionele velden

- Low and High alarms (Alarmen Laag en Hoog): Voer de BPCdrukwaarden in waarmee de alarmen High (Hoog) en Low (Laag) worden geactiveerd. Deze alarmen hebben geen invloed op lopende processen, maar waarschuwen gebruikers dat bepaalde omstandigheden buiten de gespecificeerde gebruiksparameters vallen.
- Low Low and High High alarms (Alarmen Laag-laag en Hooghoog): Vul de velden voor de interlockvergelijking en hysterese in om de voorwaarden in te stellen waarmee de alarmen voor BPC-druk High High (Hoog-hoog) en Low Low (Laag-laag) worden geactiveerd. Bij activering van deze alarmen klinkt een zoemer en worden lopende processen gepauzeerd of gestopt, afhankelijk van de gebruikersinstellingen. Voor de alarmen High High (Hoog-hoog) en Low Low (Laag-laag) moeten gebruikers waarden opgeven in de velden 'Interlock' en 'Hysteresis' (Hysterese). Deze waarden bepalen onder welke voorwaarden een alarm wordt geactiveerd, welke acties vervolgens moeten worden ondernomen en wanneer de functies worden hervat.
- Calbration (Kalibratie): Configureer een nieuwe kalibratie voor BPC-druk met een éénpunts offset. Zie hoofdstuk 4 voor meer informatie over kalibratieprocedures.

 Zoek de module 'BPC Pressure' (BPC-druk) op het startscherm van de touchscreenconsole op en raak 'Configure' (Configureren) aan. Het scherm BPC Module Configuration (BPC-module configureren) verschijnt (afbeelding 3.25). Raak het veld 'Units' (Eenheden) aan.

BPC Module Configuration				
Units:	PSI	~		
Low and High Alarms:	Lo Disabled-Hi Disab	•		
Low Low and High High Alarms:	LoLo Disabled-HiHi	•		
Alarm Designation (Pumps):	LoLo- None, HiHi- N	•		
Calibration:	Calibrate	▼		
		Done		

Afbeelding 3.25. Scherm BPC Module Configuration (BPC-module configureren).

- Selecteer mbar of psi in het geopende scherm. Raak 'Next' (Volgende) aan. Het scherm BPC Pressure Configuration (BPCdruk configureren) wordt opnieuw geopend. De geselecteerde eenheden worden weergegeven in het veld naast 'Units' (Eenheden). Raak vervolgens het veld 'Maximum air pressure value' (Maximale luchtdrukwaarde) aan.
- 3. Als u de alarmen High (Hoog)/Low (Laag) wilt instellen voor luchtdruk, raakt u het veld 'Low and High alarms' (Alarmen Laag en Hoog) aan. Het geopende scherm bevat aparte gedeeltes voor de alarmen High (Hoog) en Low (Laag) (afbeelding 3.26). Als u het alarm High (Hoog) wilt instellen, schuift u de knop On/Off (Aan/ Uit) naar de kant 'On' (Aan). Voer in het veld bij 'High Alarm' (Hoog alarm) de luchtdrukwaarde in waarbij het alarm High (Hoog) wordt geactiveerd. Raak 'Next' (Volgende) aan.

۲	BPC Pressure High and Low Alarms		
	Would you like to set the H	igh and Low alar	ms?
		High Alarm	
	Off On	0.00	
		PSI	
		Low Alarm	
	Off On On	0.00	
		PSI	
		Cancel	Done

Afbeelding 3.26. Scherm BPC Pressure High and Low Alarms (Alarmen Hoog en Laag voor BPC-druk).

- Als u het alarm Low (Laag) wilt instellen, schuift u de knop On/Off (Aan/Uit) naar de kant 'On' (Aan). Voer in het veld bij 'Low Alarm' (Laag alarm) de luchtdrukwaarde in waarbij het alarm Low (Laag) wordt geactiveerd.
- Nadat u de waarden voor de alarmen High (Hoog) en Low (Laag) hebt ingevoerd, raakt u 'Done' (Gereed) aan. In het scherm BPC Pressure Configuration (BPC-druk configureren) worden nu de ingevoerde waarden voor de alarmen High (Hoog) en Low (Laag) weergegeven.
- Als u de alarmen High High (Hoog-hoog)/Low Low (Laag-laag) wilt instellen voor luchtdruk in verband met pompen, raakt u het veld 'Low Low and High High alarms' (Alarmen Laag-laag en Hooghoog) aan.
- 7. Voor de alarmen High High (Hoog-hoog) en Low Low (Laaglaag) worden er velden met het label 'Interlock' en 'Hysteresis' (Hysterese) weergegeven. Raak 'Interlock' in een van de alarmvelden aan om de initiële voorwaarde in te stellen waarbij het alarm wordt geactiveerd, of 'Hysteresis' (Hysterese) om aan te geven welke actie na activering moet worden uitgevoerd.
- In het geopende scherm worden de interlock- en hysteresegedeeltes getoond (afbeelding 3.27). In het gedeelte 'Agitation Interlock Equation' (Interlockvergelijking voor roeren) voert u waarden in bij Control Set Point (Controle-instelpunt) (in mbar of psi) en selecteert u een waarde bij Action Taken (Ondernomen actie) (Pause (Pauzeren) of Stop (Stoppen)).



Afbeelding 3.27. Scherm BPC Pressure High High and Low Low Alarms (Alarmen Hoog-hoog en Laag-laag voor BPC-druk).

- 9. Het gedeelte 'Hysteresis' (Hysterese) (onder het gedeelte Interlock) bepaalt de luchtdruk die de BPC moet bereiken (en hoe lang die luchtdruk moet worden aangehouden, in seconden) voordat processen worden hervat. Voer een waarde in bij Set Point (Instelpunt) (in mbar of psi) en Seconds (Seconden). Wanneer u gereed bent, raakt u 'Done' (Gereed) aan.
- 10. Nadat u de interlock- en hystereseparameters voor de alarmen High High (Hoog-hoog) en Low Low (Laag-laag) hebt ingesteld, raakt u 'Done' (Gereed) aan. In het opnieuw geopende scherm BPC Pressure Configuration (BPC-druk configureren) worden nu de ingevoerde waarden voor de alarmen High High (Hoog-hoog) en Low Low (Laag-laag) weergegeven.
- 11. Als u een kalibratie voor BPC-druk wilt instellen, raakt u het veld 'Calibration' (Kalibratie) aan. Volg de kalibratie-instructies uit hoofdstuk 4 — Kalibratieprocedures. U kunt de kalibratie ook overslaan en 'Done' (Gereed) aanraken om terug te gaan naar het startscherm.

3.5.6 De module Pumps (Pompen) configureren

Gebruik de volgende informatie en stappen om de module Pumps (Pompen) te configureren. Elk van de onderstaande velden wordt geïllustreerd in afbeelding 3.28. **Opmerking:** U kunt 'Done' (Gereed) onder aan het scherm Configuration (Configuratie) te allen tijde aanraken tijdens de configuratie van de module. Uw invoer wordt opgeslagen en u kunt doorgaan waar u was gebleven wanneer u de moduleconfiguratie opnieuw opent.

Verplichte velden

- **RPM scaling (RPM-schaling):** Zet het uitvoersignaal voor de pomp om van milliampères (mA) in RPM.
- **RPM limits (RPM-limieten):** Stel de maximale en minimale RPMwaarden voor de pomp in.

Optionele velden

- Pump name (Naam van pomp): Elke pompmodule (maximaal vier) kan een aparte naam krijgen om het doel van de pompen aan te geven. Eén pomp zou bijvoorbeeld de naam 'P1 zuur' kunnen krijgen en een andere de naam 'P2 aftappen' zodat gebruikers de pompen gemakkelijker van elkaar kunt onderscheiden.
 Opmerking: Alle pompnamen krijgen een voorvoegsel 'P#' om aan te geven welke pomppoort op de pomp aangesloten is.
- **Calibration (Kalibratie):** Configureer een nieuwe pomp met vierpunts kalibratie en met massa- of volume-eenheden. Zie hoofdstuk 4 voor meer informatie over kalibratieprocedures.
- Zoek de module Pump (Pomp) op het startscherm van de touchscreenconsole op en raak 'Configure' (Configureren) aan. Het scherm Pump Configuration (Pomp configureren) verschijnt (afbeelding 3.28). Raak het veld 'RPM Scaling' (RPM-schaling) aan.

Pump1 Module Configuration				
Module Name:	Pump1	~		
RPM Scaling:	4.00mA=0 20.00mA=			
RPM Limits:	0-300	•		
Calibration:	Not Calibrated	~		
		Done		

Afbeelding 3.28. Scherm Pump Configuration (Pomp configureren).

 In het geopende scherm (afbeelding 3.29) wordt u gevraagd de RPM-waarden op te geven die u op de touchscreenconsole wilt weergeven na omzetting van het uitvoersignaal (in mA). Raak 'Done' (Gereed) aan wanneer u klaar bent.

Pump1 RPM Scaling				
	Convert Outpu 20.00 Max. mA	it Sigr = [nal for Pump1 300 Max. RPM	
	4.00 Min. mA		0 Min. RPM	
			Cancel	Done

Afbeelding 3.29. Scherm Pump RPM Scaling (RPM-schaling pomp).

- Het scherm Pump Configuration (Pomp configureren) wordt opnieuw geopend. De ingevoerde schalingswaarden worden weergegeven in het veld naast 'RPM scaling' (RPM-schaling). Raak vervolgens het veld 'RPM Limits' (RPM-limieten) aan.
- Stel in het geopende scherm de RPM-limieten van de pomp in door de maximum- en minimumsnelheden (in RPM) op te geven. Raak 'Done' (Gereed) aan.
- 5. Als u de pompmodule een naam wilt geven raakt u het veld 'Pump name' (Naam van pomp) aan. Geef een naam op voor de pomp met het alfanumerieke toetsenbord dat op het scherm verschijnt. Wanneer u gereed bent, raakt u 'Enter' (Invoeren) aan. Het scherm Pump Configuration (Pomp configureren) wordt opnieuw geopend met de naam die u naast het veld 'Pump Name' (Naam van pomp) hebt opgegeven.
- Als u de pomp wilt kalibreren, raakt u het veld 'Calibration' (Kalibratie) aan. Volg de kalibratie-instructies uit hoofdstuk 4 — Kalibratieprocedures. U kunt de kalibratie ook overslaan en 'Done' (Gereed) aanraken om terug te gaan naar het startscherm.

3.5.7 De module Fill (Vullen) configureren

Gebruik de volgende informatie en stappen om de module Fill (Vullen) te configureren. Elk van de onderstaande velden wordt geïllustreerd in afbeelding 3.30. De module Mass (Massa) moet zijn geconfigureerd voordat u de module Fill (Vullen) configureert. **Opmerking:** U kunt 'Done' (Gereed) onder aan het scherm Configuration (Configuratie) te allen tijde aanraken tijdens de configuratie van de module. Uw invoer wordt opgeslagen en u kunt doorgaan waar u was gebleven wanneer u de moduleconfiguratie opnieuw opent.

Verplichte velden

- **Pinch valve (Knijpventiel):** Dit veld wordt weergegeven als de touchscreenconsole wel of geen knijpventiel heeft gedetecteerd.
- Units and liquid density (Eenheden en vloeistofdichtheid): Selecteer de eenheden die u wilt gebruiken (liter (l) of kilogram (kg)).
- Fill pump selection (Vulpomp selecteren): Selecteer een pomp om aan de module Fill (Vullen) toe te wijzen. Als u niet onmiddellijk een pomp wilt toewijzen, raakt u 'Assign Pump Later' (Pomp later toewijzen) aan.
- Fill control (Vullen reguleren): Stel de parameters in voor het vulproces, met inbegrip van de totale vloeistoftoediening, stroomsnelheden faseren, overgang tussen fasen en beëindiging.

Optionele velden

- Geen
- Zoek de module Fill (Vullen) op het startscherm van de touchscreenconsole op en raak 'Configure' (Configureren) aan. Het scherm Fill Module Configuration (Vulmodule configureren) verschijnt (afbeelding 3.30). Raak het veld 'Units and Liquid Density' (Eenheden en vloeistofdichtheid) aan.

\odot	Fill Module Configuration		
	Fill Pinch valve:	Detected	
	Units and Liquid Density:	kg, 1.000 g/mL	•
	Fill Pump Selection:		•
	Fill Control:	Setpoint 25.0 kg	•
			Done

Afbeelding 3.30. Scherm Fill Module Configuration (Vulmodule configureren).

 Selecteer in het geopende scherm de eenheden die u wilt gebruiken (Mass (Massa) (in kg) of Volume (in l)). Geef vervolgens de vloeistofdichtheid op (in g/ml). Raak 'Done' (Gereed) aan wanneer u klaar bent.

- Het scherm Fill Module Configuration (Vulmodule configureren) wordt opnieuw geopend. De geselecteerde eenheden en vloeistofdichtheid worden weergegeven in het veld naast 'Units and Liquid Density' (Eenheden en vloeistofdichtheid). Raak vervolgens het veld 'Fill pump selection' (Vulpomp selecteren) aan.
- 4. In het geopende scherm wordt u gevraagd een pomp te selecteren voor toewijzing aan de module Fill (Vullen). Nadat u de pomp hebt aangeraakt die u voor het vullen wilt gebruiken, raakt u 'Done' (Gereed) aan. **Opmerking:** Er moet ten minste één pomp op het systeem zijn aangesloten om een pomp aan de module Fill (Vullen) toe te wijzen.
- Het scherm Fill Configuration (Vulmodule configureren) wordt opnieuw geopend, met de pomp die u naast het veld 'Fill Pump Selection' (Vulpomp selecteren) hebt geselecteerd. Raak vervolgens 'Fill control' (Vullen reguleren) aan.
- 6. Het geopende scherm bevat velden voor de verschillende stappen van het vulproces. Voer de gewenste waarden in voor de volgende velden:
 - **Total fluid delivery (Totale vloeistoftoediening):** De totale hoeveelheid vloeistof die u in de BPC wilt pompen.
 - Stage 1 flow rate (Flowsnelheid fase 1): De snelheid waarmee u de BPC in eerste instantie wilt vullen (in I of kg per minuut).
 - Stage 1 to Stage 2 transition (Overgang fase 1 naar fase 2): Het percentage vloeistof dat in de BPC moet worden gepompt voordat de pomp naar fase 2 gaat.
 - Stage 2 flow rate (Flowsnelheid fase 2): De gewenste snelheid bij voltooiing van het vulproces. De stroomsnelheid van fase 2 moet aanzienlijk lager zijn dan die van fase 1 om overlopen te voorkomen.
 - Stage 2 termination (Beëindiging fase 2): De actie die de pompen moeten uitvoeren wanneer het vulproces voltooid is (Pause (Pauzeren) of Stop (Stoppen)).
- 7. Raak 'Done' (Gereed) aan om terug te gaan naar het startscherm.

3.5.8 De module Harvest (Aftappen) configureren

Gebruik de volgende informatie en stappen om de module Harvest (Oogsten) te configureren. Elk van de onderstaande velden wordt geïllustreerd in afbeelding 3.31. De module Mass (Massa) moet zijn geconfigureerd voordat u de module Harvest (Aftappen) configureert. **Opmerking:** U kunt 'Done' (Gereed) onder aan het scherm Configuration (Configuratie) te allen tijde aanraken tijdens de configuratie van de module. Uw invoer wordt opgeslagen en u kunt doorgaan waar u was gebleven wanneer u de moduleconfiguratie opnieuw opent.

Verplichte velden

- **Pinch valve (Knijpventiel):** Dit veld wordt weergegeven als de touchscreenconsole wel of geen knijpventiel heeft gedetecteerd.
- Units and liquid density (Eenheden en vloeistofdichtheid): Selecteer de eenheden die u wilt gebruiken (liter (l) of kilogram (kg)).
- Harvest pump selection (Aftappomp selecteren): Selecteer een pomp om aan de module Harvest (Aftappen) toe te wijzen. Als u niet onmiddellijk een pomp wilt toewijzen, raakt u 'Assign Pump Later' (Pomp later toewijzen) aan.
- Harvest control (Aftappen reguleren): Stel de parameters in voor het aftapproces, met inbegrip van de totale vloeistoftoediening, stroomsnelheden faseren, overgang tussen fasen en beëindiging.

Optionele velden

- Geen
- Zoek de module Harvest (Aftappen) op het startscherm van de touchscreenconsole op en raak 'Configure' (Configureren) aan. Het scherm Harvest Module Configuration (Aftapmodule configureren) verschijnt (afbeelding 3.31). Raak het veld 'Units and Liquid Density' (Eenheden en vloeistofdichtheid) aan.

lacksquare	Harvest Module Configuration		
	Harvest Pinch valve:	Detected	
	Units and Liquid Density:	kg, 1.000 g/mL	~
	Harvest Pump Selection:		~
	Harvest Control:	Setpoint 0.0 kg	~
			Done

Afbeelding 3.31. Scherm Harvest Module Configuration (Aftapmodule configureren).

 Selecteer in het geopende scherm de eenheden die u wilt gebruiken (Mass (Massa) (in kg) of Volume (in l)). Geef vervolgens de vloeistofdichtheid op (in g/ml). Raak 'Done' (Gereed) aan wanneer u klaar bent.

- Het scherm Harvest Module Configuration (Aftapmodule configureren) wordt opnieuw geopend. De geselecteerde eenheden en vloeistofdichtheid worden weergegeven in het veld naast 'Units and Liquid Density' (Eenheden en vloeistofdichtheid). Raak vervolgens het veld 'Harvest pump selection' (Aftappomp selecteren) aan.
- 4. In het geopende scherm wordt u gevraagd een pomp te selecteren voor toewijzing aan de module Harvest (Aftappen). Nadat u de pomp hebt aangeraakt die u voor het aftappen wilt gebruiken, raakt u 'Done' (Gereed) aan. **Opmerking:** Er moet ten minste één pomp op het systeem zijn aangesloten om een pomp aan de module Harvest (Aftappen) toe te wijzen.
- 5. Het scherm Harvest Module Configuration (Aftapmodule configureren) wordt opnieuw geopend , met de pomp die u naast het veld 'Harvest Pump Selection' (Aftappomp selecteren) hebt geselecteerd. Raak vervolgens 'Harvest Control' (Aftappen reguleren) aan.
- 6. Het geopende scherm bevat velden voor de verschillende stappen van het aftapproces. Voer de gewenste waarden in voor de volgende velden:
 - **Total fluid delivery (Totale vloeistoftoediening):** De totale hoeveelheid vloeistof die u uit de BPC wilt pompen.
 - Stage 1 flow rate (Flowsnelheid fase 1): De snelheid waarmee u de BPC in eerste instantie wilt aftappen (in I of kg per minuut).
 - Stage 1 to Stage 2 transition (Overgang fase 1 naar fase 2): Het percentage vloeistof dat uit de BPC moet worden gepompt voordat de pomp naar fase 2 gaat.
 - Stage 2 flow rate (Flowsnelheid fase 2): De gewenste snelheid bij voltooiing van het aftapproces. De stroomsnelheid van fase 2 moet aanzienlijk lager zijn dan die van fase 1 om overmatig aftappen te voorkomen.
 - Stage 2 termination (Beëindiging fase 2): De actie die de pompen moeten uitvoeren wanneer het aftapproces voltooid is (Pause (Pauzeren) of Stop (Stoppen)).
- 7. Raak 'Done' (Gereed) aan om terug te gaan naar het startscherm.

3.5.9 De pH-module configureren

Gebruik de volgende informatie en stappen om de pH-module te configureren. Elk van de onderstaande velden wordt geïllustreerd in afbeelding 3.32. **Opmerking:** U kunt 'Done' (Gereed) onder aan het scherm Configuration (Configuratie) te allen tijde aanraken tijdens de configuratie van de module. Uw invoer wordt opgeslagen en u kunt doorgaan waar u was gebleven wanneer u de moduleconfiguratie opnieuw opent.

Verplichte velden

- pH sensors available (Beschikbare pH-sensoren): Selecteer welke pH-sensoren u voor gebruik beschikbaar wilt maken. U kunt pH1, pH2 of beide selecteren.
- Active pH sensor (Actieve pH-sensor): Selecteer welke pHsensor u actief wilt maken. U kunt pH1 of pH2 selecteren.

Optionele velden

- Temperature compensation (Temperatuurcompensatie): Als de module Temperature (Temperatuur) is geconfigureerd, wordt de aflezing van de temperatuursensor gebruikt voor pH-temperatuurcompensatie. Als de module Temperature (Temperatuur) niet is geconfigureerd, voert u de Expected Vessel Temperature (Verwachte vattemperatuur) en wordt deze gebruikt. U kunt er ook voor kiezen de module Temperature (Temperatuur) te configureren voordat u doorgaat met de pH-configuratie.
- Low and High alarms (Alarmen Laag en Hoog): Voer de pHwaarden in waarmee de pH-alarmen High (Hoog) en Low (Laag) worden geactiveerd. Deze alarmen hebben geen invloed op lopende processen, maar waarschuwen gebruikers dat bepaalde omstandigheden buiten de gespecificeerde gebruiksparameters vallen.
- Low Low and High High alarms (Alarmen Laag-laag en Hooghoog): Vul de velden voor de interlockvergelijking en hysterese in om de voorwaarden in te stellen waarmee de pH-alarmen High High (Hoog-hoog) en Low Low (Laag-laag) worden geactiveerd. Bij activering van deze alarmen klinkt een zoemer en worden lopende processen gepauzeerd of gestopt, afhankelijk van de gebruikersinstellingen. Voor de alarmen High High (Hoog-hoog) en Low Low (Laag-laag) moeten gebruikers waarden opgeven in de velden 'Interlock' en 'Hysteresis' (Hysterese). Deze waarden bepalen onder welke voorwaarden een alarm wordt geactiveerd, welke acties vervolgens moeten worden ondernomen en wanneer de functies worden hervat.
- Alarm designation (Alarmaanduiding): Nadat u de pH-alarmen High High (Hoog-hoog) en/of Low Low (Laag-laag) hebt ingesteld, gebruikt u dit veld om te selecteren op welke pompen u deze alarmen wilt toepassen. Bijvoorbeeld voor Pomp1 (Zuur) schakelt u het alarm High High (Hoog-hoog) in en het alarm Low Low (Laag-laag) uit.
- Calibration (Kalibratie): Configureer een nieuwe pH-kalibratie met directe invoer, 2-punts of 3-punts kalibratie. Zie hoofdstuk 4 voor meer informatie over kalibratieprocedures.
- Zoek de pH-module op het startscherm van de touchscreenconsole op en raak 'Configure' (Configureren) aan. Het scherm pH Module Configuration (pH-module configureren) verschijnt (afbeelding 3.32). Raak het veld 'pH sensors available' (Beschikbare pH-sensoren) aan

PH Module Configuration			
nH Sensors Available:	pH1 and pH2		
Active pH Sensor:	pH1	~	
Temperature Compensation:	Temp1 (45.00°C)	•	
Low and High Alarms:	Lo Disabled-Hi Disab	•	
Low Low and High High Alarms:	LoLo Disabled-HiHi	•	
Alarm Designation (Pumps):	LoLo- None, HiHi- N	•	
		Done	

Afbeelding 3.32. Scherm pH Module Configuration (pH-module configureren).

- 2. Selecteer in het geopende scherm de pH-sensor(en) die u voor gebruik beschikbaar wilt maken. U kunt pH1, pH2 of beide selecteren. Raak 'Done' (Gereed) aan wanneer u klaar bent.
- Het scherm pH Module Configuration (pH-module configureren) wordt opnieuw geopend. De geselecteerde sensoren worden weergegeven in het veld naast 'pH Sensors Available' (Beschikbare pH-sensoren). Raak vervolgens het veld 'Active pH sensor' (Actieve pH-sensor) aan.

- Selecteer in het geopende scherm de pH-sensor die u wilt activeren. U kunt pH1 of pH2 selecteren. Wanneer u de actieve sensor hebt geselecteerd, raakt u 'Done' (Gereed) aan.
 Opmerking: U kunt te allen tijde een andere sensor activeren door terug te gaan naar het scherm pH Module Configuration (pH-module activeren) en Active pH sensor (Actieve pH-sensor) te selecteren.
- 5. Het scherm pH Module Configuration (pH-module configureren) wordt opnieuw geopend. Raak het veld 'Temperature Compensation' (Temperatuurcompensatie) aan.
 - Als de module Temperature (Temperatuur) niet is geconfigureerd, wordt u in een scherm gevraagd de module Temperature (Temperatuur) te configureren (om de aflezing van de temperatuursensor te gebruiken) of de verwachte vattemperatuur handmatig op te geven.
 - Als de module Temperature (Temperatuur) is geconfigureerd, kunt u kiezen de temperatuursensor in of uit te schakelen. Als u ervoor kiest de sensor uit te schakelen, moet u de verwachte vattemperatuur handmatig opgeven (afbeelding 3.33).

	pH Temperature Compensation
т	he Expected Vessel Temperature will be used if there is no RTD present
	Expected Vessel Temperature
	25.00
	°C
	You can use the RTD if present by setting the slider to enable below
	OffO On 45.00 °C
	Cancel Done

Afbeelding 3.33. Scherm pH Temperature Compensation (pH-temperatuurcompensatie).

Wanneer u de module Temperature (Temperatuur) hebt geconfigureerd en/of hebt ingeschakeld of de verwachte vattemperatuur hebt opgegeven, raakt u 'Done' (Gereed) aan.

- Het scherm pH Module Configuration (pH-module configureren) wordt opnieuw geopend. Als u de alarmen High (Hoog)/Low (Laag) wilt instellen voor pH, raakt u het veld 'Low and High alarms' (Alarmen Laag en Hoog) aan.
- Het geopende scherm bevat aparte gedeeltes voor de alarmen High (Hoog) en Low (Laag). Als u het alarm High (Hoog) wilt configureren, schuift u de knop On/Off (Aan/Uit) naar de kant 'On' (Aan). Voer in het veld bij 'High Alarm' (Hoog alarm) de pHwaarde in waarbij het alarm High (Hoog) wordt geactiveerd.
- Als u het alarm Low (Laag) wilt configureren, schuift u de knop On/Off (Aan/Off) naar de kant 'On' (Aan). Voer in het veld bij 'Low Alarm' (Laag alarm) de pH-waarde in waarbij het alarm Low (Laag) wordt geactiveerd.
- 9. Nadat u de waarden voor de alarmen High (Hoog) en Low (Laag) hebt ingevoerd, raakt u 'Done' (Gereed) aan.
- 10. In het scherm pH Configuration (pH configureren) worden nu de ingevoerde waarden voor de alarmen High (Hoog) en Low (Laag) weergegeven. Als u de alarmen High High (Hoog-hoog)/Low Low (Laag-laag) wilt instellen voor pH, raakt u het veld 'Low Low and High High alarms' (Alarmen Laag-laag en Hoog-hoog) aan.
- Het geopende scherm (afbeelding 3.34) bevat aparte gedeeltes voor de alarmen High High (Hoog-hoog) en Low Low (Laag-laag). Schuif de knoppen On/Off (Aan/uit) in elk gedeelte naar de kant 'On' (Aan) om de alarmen in te schakelen.



Afbeelding 3.34. Scherm High High and Low Low Alarms (Alarmen Hoog-hoog en Laag-laag).

- 12. Voor de alarmen High High (Hoog-hoog) en Low Low (Laaglaag) worden er velden met het label 'Interlock' en 'Hysteresis' (Hysterese) weergegeven. Raak 'Interlock' in een van de alarmvelden aan om de initiële voorwaarde in te stellen waarbij het alarm wordt geactiveerd, en 'Hysteresis' (Hysterese) om aan te geven welke actie na activering moet worden uitgevoerd.
- In het geopende scherm worden de interlock- en hysteresegedeeltes getoond. In het gedeelte 'Interlock Equation' (Interlockvergelijking) voert waarden in bij Control Set Point (Controle-instelpunt) (in pH) en selecteert u een waarde bij Action Taken (Ondernomen actie) (Pause (Pauzeren) of Stop (Stoppen)).
- 14. Het gedeelte 'Hysteresis' (Hysterese) (onder het gedeelte Interlock) bepaalt wanneer processen worden hervat nadat het alarm is geactiveerd. Voer de tolerantie en de tijdsduur in gedurende welke het pH-niveau van het controle-instelpunt (plus de tolerantie) moet worden vastgehouden voordat wordt verdergegaan (in seconden). Wanneer u gereed bent, raakt u 'Done' (Gereed) aan.
- 15. Wanneer u de interlock en hysterese voor het alarm High High (Hoog-hoog) en Low Low (Laag-laag) hebt ingesteld, raakt u 'Done' (Gereed) aan in het scherm 'pH High High and Low Low Alarms' (Alarmen Hoog-hoog en Laag-laag voor pH).
- 16. Als u alarmen voor specifieke, voor de S.U.M. geconfigureerde pompen wilt aanduiden, raakt u het veld 'Alarm Designation' (Alarmaanduiding) in het scherm pH Configuration (pHconfiguratie) aan. **Opmerking:** Wanneer een pomp is ingeschakeld, kan het pH-niveau worden aangepast door na activering van het alarm een bolus toe te dienen.
- 17. In het geopende scherm (afbeelding 3.35) ziet u de interlock- en hystereseparameters die u hebt opgegeven voor de alarmen High High (Hoog-hoog) en Low Low (Laag-laag). Eronder worden de met het systeem ingestelde pompen getoond. Raak een pomp aan om de alarmen High High (Hoog-hoog) en Low Low (Laaglaag) in of uit te schakelen.



Afbeelding 3.35. Scherm pH Pump Designation (pH-pompaanduiding).

- 18. Schuif op het geopende scherm de knop Enable/Disable (In-/ uitschakelen) naar de kant Enable (Inschakelen) voor de alarmen High High (Hoog-hoog) en/of Low Low (Laag-laag) om het of de alarmen in te schakelen voor de geselecteerde pomp. Herhaal deze stap om alarmen toe te wijzen aan alle andere geconfigureerde pompen. Wanneer u gereed bent, raakt u de knop 'Done' (Gereed) aan.
- 19. Als u de pH wilt kalibreren, raakt u het veld 'Calibration' (Kalibratie) aan. Volg de kalibratie-instructies uit hoofdstuk 4 — Kalibratieprocedures. U kunt de kalibratie ook overslaan en 'Done' (Gereed) aanraken om terug te gaan naar het startscherm.

3.5.10 De module Conductivity (Geleidbaarheid) configureren

Gebruik de volgende informatie en stappen om de module Conductivity (Geleidbaarheid) te configureren. Elk van de onderstaande velden wordt geïllustreerd in afbeelding 3.36. **Opmerking:** U kunt 'Done' (Gereed) onder aan het scherm Configuration (Configuratie) te allen tijde aanraken tijdens de configuratie van de module. Uw invoer wordt opgeslagen en u kunt doorgaan waar u was gebleven wanneer u de moduleconfiguratie opnieuw opent.

Verplichte velden

- Units (Eenheden): Selecteer de eenheden die u wilt gebruiken (μS/cm of mS/cm).
- Active conductivity sensor (Actieve geleidbaarheidssensor): Selecteer welke geleidbaarheidssensor u actief wilt maken. U kunt Conductivity1 of Conductivity2 selecteren.

Optionele velden

- Low and High alarms (Alarmen Laag en Hoog): Voer de geleidbaarheidswaarden in waarmee de geleidbaarheidsalarmen High (Hoog) en Low (Laag) worden geactiveerd. Deze alarmen hebben geen invloed op lopende processen, maar waarschuwen gebruikers dat bepaalde omstandigheden buiten de gespecificeerde gebruiksparameters vallen.
- Low Low and High High alarms (Alarmen Laag-laag en Hoog-hoog): Vul de velden voor de interlockvergelijking en hysterese in om de voorwaarden in te stellen waarmee de geleidbaarheidsalarmen High High (Hoog-hoog) en Low Low (Laag-laag) worden geactiveerd. Bij activering van deze alarmen klinkt een zoemer en worden lopende processen gepauzeerd of gestopt, afhankelijk van de gebruikersinstellingen. Voor de alarmen High High (Hoog-hoog) en Low Low (Laag-laag) moeten gebruikers waarden opgeven in de velden 'Interlock' en 'Hysteresis' (Hysterese). Deze waarden bepalen onder welke voorwaarden een alarm wordt geactiveerd, welke acties vervolgens moeten worden ondernomen en wanneer de functies worden hervat.
- Alarm designation (Alarmaanduiding): Nadat u de geleidbaarheidsalarmen High High (Hoog-hoog) en/of Low Low (Laag-laag) hebt ingesteld, gebruikt u dit veld om te selecteren op welke pompen u deze alarmen wilt toepassen. Bijvoorbeeld voor Pomp1 schakelt u het alarm Low Low (Laag-laag) in en het alarm High High (Hoog-hoog) uit.
- Calibration (Kalibratie): Configureer een nieuwe geleidbaarheidskalibratie met directe invoer, 2-punts of 3-punts kalibratie. Zie hoofdstuk 4 voor meer informatie over kalibratieprocedures.
- Zoek de module Conductivity (Geleidbaarheid) op het startscherm van de touchscreenconsole op en raak 'Configure' (Configureren) aan. Het scherm Conductivity Module Configuration (Module Geleidbaarheid configureren) verschijnt (afbeelding 3.36). Raak het veld 'Units' (Eenheden) aan.

Conductivity Module Configuration			
Units:	µS/cm	•	
Active Conductivity Sensor:	Conductivity1	•	
Low and High Alarms:	Lo Disabled-Hi Disab	•	
Low Low and High High Alarms:	LoLo Disabled-HiHi	•	
Alarm Designation (Pumps):	LoLo- None, HiHi- N	•	
Calibration:	Calibrate	•	
		Done	

Afbeelding 3.36. Scherm Conductivity Configuration (Geleidbaarheid configureren).

- Selecteer in het geopende scherm de eenheden die u voor de touchscreenconsole beschikbaar wilt maken. U kunt μS/cm of mS/cm selecteren. Raak 'Next' (Volgende) aan.
- Het scherm Conductivity Module Configuration (Module Geleidbaarheid configureren) wordt opnieuw geopend. Raak 'Active Conductivity Sensor' (Actieve geleidbaarheidssensor) aan. In het geopende scherm wordt gevraagd welke geleidbaarheidssensor de gebruiker wil activeren. U kunt Conductivity1 of Conductivity2 selecteren. Raak 'Next' (Volgende) aan.
- Het scherm Conductivity Module Configuration (Module Geleidbaarheid configureren) wordt opnieuw geopend. Als u de alarmen High (Hoog)/Low (Laag) wilt instellen voor geleidbaarheid, raakt u het veld 'Low and High Alarms' (Alarmen Laag en Hoog) aan.
- Het geopende scherm bevat aparte gedeeltes voor de alarmen High (Hoog) en Low (Laag) (afbeelding 3.37). Als u het alarm High (Hoog) wilt configureren, schuift u de knop On/Off (Aan/ Uit) naar de kant 'On' (Aan). Voer in het veld 'High Alarm' (Hoog alarm) het geleidbaarheidsniveau in waarbij het alarm High (Hoog) wordt geactiveerd.

۲	Co	nductivity High a	nd Low Alarms	
Would you like to set the High and Low alarms?				
			High Alarm	
	Off	On	0	
			µS/cm	
			Low Alarm	
	Off	On On	0	
			µS/cm	
			Cancel	Done

Afbeelding 3.37. Conductivity High and Low Alarms (Alarmen Hoog en Laag voor geleidbaarheid).

- Als u het alarm Low (Laag) wilt configureren, schuift u de knop On/Off (Aan/Off) naar de kant 'On' (Aan). Voer in het veld 'Low Alarm' (Laag alarm) het geleidbaarheidsniveau in waarbij het alarm Low (Laag) wordt geactiveerd.
- 7. Nadat u de waarden voor de alarmen High (Hoog) en Low (Laag) hebt ingevoerd, raakt u 'Done' (Gereed) aan.
- 8. In het scherm Conductivity Module Configuration (Module Geleidbaarheid configureren) worden nu de ingevoerde waarden voor de alarmen High (Hoog) en Low (Laag) weergegeven. Als u de alarmen High High (Hoog-hoog)/Low Low (Laag-laag) wilt instellen voor, raakt u het veld 'Low Low and High High alarms' (Alarmen Laag-laag en Hoog-hoog) aan.
- Het geopende scherm (afbeelding 3.38) bevat aparte gedeeltes voor de alarmen High High (Hoog-hoog) en Low Low (Laag-laag). Schuif de knoppen On/Off (Aan/uit) in elk gedeelte naar de kant 'On' (Aan) om de alarmen in te schakelen.

€	Conductivity1 High High and Low Low Alarms		
	High High Alarm	Off O	
	Interlock:	Conductivity1>0µS/cm then Stop	
	Hysteresis:	Resume if Conductivity1<0µS/cm for 0s	
	Low Low Alarm	Off O	
	Interlock:	Conductivity1<0µS/cm then Stop	
	Hysteresis:	Resume if Conductivity1>0µS/cm for 0s	
		Cancel Done	

Afbeelding 3.38. Scherm High High and Low Low Alarms (Alarmen Hoog-hoog en Laag-laag)

- 10. Voor de alarmen High High (Hoog-hoog) en Low Low (Laaglaag) worden er velden met het label 'Interlock' en 'Hysteresis' (Hysterese) weergegeven. Raak 'Interlock' in een van de alarmvelden aan om de initiële voorwaarde in te stellen waarbij het alarm wordt geactiveerd, en 'Hysteresis' (Hysterese) om aan te geven welke actie na activering moet worden uitgevoerd.
- 11. In het geopende scherm worden de interlock- en hysteresegedeeltes getoond. In het gedeelte 'Interlock Equation' (Interlockvergelijking) voert u waarden in bij Control Set Point (Controle-instelpunt) en selecteert u een waarde bij Action Taken (Ondernomen actie) (Pause (Pauzeren) of Stop (Stoppen)).
- 12. Het gedeelte 'Hysteresis' (Hysterese) (onder het gedeelte Interlock) bepaalt wanneer processen worden hervat nadat het alarm is geactiveerd. Voer de tolerantie en de tijdsduur in gedurende welke het geleidbaarheidsniveau van het controle-instelpunt voor (plus de tolerantie) moet worden vastgehouden voordat wordt verdergegaan (in seconden). Wanneer u gereed bent, raakt u 'Done' (Gereed) aan.
- 13. Nadat u de interlock- en hystereseparameters voor de alarmen High High (Hoog-hoog) en Low Low (Laag-laag) hebt ingesteld, raakt u 'Done' (Gereed) aan.
- 14. Als u alarmen voor specifieke pompen wilt aanduiden, raakt u het veld 'Alarm Designation' (Alarmaanduiding) in het scherm Conductivity Module Configuration (Module Geleidbaarheid configureren) aan. **Opmerking:** Wanneer een pomp is ingeschakeld, kan het geleidbaarheidsniveau worden aangepast door na activering van het alarm een bolus toe te dienen.

15. In het geopende scherm (afbeelding 3.39) ziet u de interlocken hystereseparameters die u hebt opgegeven voor de alarmen High High (Hoog-hoog) en Low Low (Laag-laag). Eronder worden de met het systeem ingestelde pompen getoond. Raak een pomp aan om de alarmen High High (Hoog-hoog) en Low Low (Laag-laag) in of uit te schakelen.



Afbeelding 3.39. Scherm Alarm Designation (Alarmaanduiding).

- 16. Schuif op het geopende scherm de knop Enable/Disable (In-/ uitschakelen) naar de kant Enable (Inschakelen) voor de alarmen High High (Hoog-hoog) en/of Low Low (Laag-laag) om het of de alarmen in te schakelen voor de geselecteerde pomp. Herhaal deze stap om alarmen toe te wijzen aan alle andere geconfigureerde pompen. Wanneer u gereed bent, raakt u de knop 'Done' (Gereed) aan.
- 17. Als u de geleidbaarheid wilt kalibreren, raakt u het veld
 'Calibration' (Kalibratie) aan. Volg de kalibratie-instructies uit hoofdstuk 4 — Kalibratieprocedures. U kunt de kalibratie ook overslaan en 'Done' (Gereed) aanraken om terug te gaan naar het startscherm.

3.5.11 De module Temperature (Temperatuur) configureren

Gebruik de volgende informatie en stappen om de module Temperature (Temperatuur) te configureren. Elk van de onderstaande velden wordt geïllustreerd in afbeelding 3.40. **Opmerking:** U kunt 'Done' (Gereed) onder aan het scherm Configuration (Configuratie) te allen tijde aanraken tijdens de configuratie van de module. Uw invoer wordt opgeslagen en u kunt doorgaan waar u was gebleven wanneer u de moduleconfiguratie opnieuw opent.

Verplichte velden

- Available temperature sensors (Beschikbare temperatuursensoren): Selecteer welke temperatuursensoren u voor gebruik beschikbaar wilt maken. U kunt Temperature 1 (Temperatuur 1), Temperature 2 (Temperatuur 2) of beide selecteren.
- Active temperature sensor (Actieve temperatuursensor): Selecteer welke geleidbaarheidssensor u actief wilt maken. U kunt Temperature1 of Temperature2 selecteren.
- TCU setpoint limits (Instelpuntlimieten TCU) (voor systemen met een TCU): Voor de maximum- en minimumtemperaturen in die u wilt gebruiken (in°C).

Optionele velden

- Low and High alarms (Alarmen Laag en Hoog): Voer de temperatuurwaarden in waarmee de temperatuuralarmen High (Hoog) en Low (Laag) worden geactiveerd. Deze alarmen hebben geen invloed op lopende processen, maar waarschuwen gebruikers dat bepaalde omstandigheden buiten de gespecificeerde gebruiksparameters vallen.
- Low Low and High High alarms (Alarmen Laag-laag en Hooghoog): Vul de velden voor de interlockvergelijking en hysterese in om de voorwaarden in te stellen waarmee de temperatuuralarmen High High (Hoog-hoog) en Low Low (Laag-laag) worden geactiveerd. Bij activering van deze alarmen klinkt een zoemer en worden lopende processen gepauzeerd of gestopt, afhankelijk van de gebruikersinstellingen. Voor de alarmen High High (Hoog-hoog) en Low Low (Laag-laag) moeten gebruikers waarden opgeven in de velden 'Interlock' en 'Hysteresis' (Hysterese). Deze waarden bepalen onder welke voorwaarden een alarm wordt geactiveerd, welke acties vervolgens moeten worden ondernomen en wanneer de functies worden hervat.
- Zoek de module Temperature (Temperatuur) op het startscherm van de touchscreenconsole op en raak 'Configure' (Configureren) aan. Het scherm Temperature Module Configuration (Module Temperatuur configureren) verschijnt (afbeelding 3.40). Raak het veld Available Temperature Sensors' (Beschikbare temperatuursensoren) aan.

Temperature Module Configuration				
Active Temperature Sensor:	Temp1	~		
TCU Set Point Limits:	4.00°C - 105.00°C	•		
Low and High Alarms:	Lo 5.00°C-Hi Disabled	~		
Low Low and High High Alarms:	LoLo Disabled-HiHi	•		
		(A)(0)()		
		Done		

Afbeelding 3.40. Scherm Temperature Module Configuration (Module Temperatuur configureren).

- Selecteer in het geopende scherm de temperatuursensor(en) die u voor gebruik beschikbaar wilt maken. U kunt Temperature1, Temperature2 of beide selecteren. Raak 'Done' (Gereed) aan wanneer u klaar bent.
- 3. Het scherm Temperature Module Configuration (Module Temperatuur configureren) wordt opnieuw geopend. De geselecteerde sensoren worden weergegeven in het veld naast 'Available Temperature Sensors' (Beschikbare temperatuursensoren). Raak vervolgens het veld 'Active Temperature Sensor' (Actieve temperatuursensor) aan.
- 4. Selecteer in het geopende scherm de temperatuursensor die u wilt activeren. U kunt Temperature1 of Temperature2 selecteren. Wanneer u de actieve sensor hebt geselecteerd, raakt u 'Done' (Gereed) aan. **Opmerking:** U kunt te allen tijde een andere sensor activeren door terug te gaan naar het scherm Temperature Module Configuration (Module Temperatuur configureren) en Active Temperature Sensor (Actieve temperatuursensor) te selecteren.
- Het scherm Temperature Module Configuration (Module Temperatuur configureren) wordt opnieuw geopend. Als u de alarmen High (Hoog)/Low (Laag) wilt instellen voor temperatuur, raakt u het veld 'Low and High alarms' (Alarmen Laag en Hoog) aan.

 Het geopende scherm (afbeelding 3.41) bevat aparte gedeeltes voor de alarmen High (Hoog) en Low (Laag). Als u het alarm High (Hoog) wilt configureren, schuift u de knop On/Off (Aan/Uit) naar de kant 'On' (Aan). Voer in het veld bij 'High Alarm' (Hoog alarm) de temperatuurwaarde in waarbij het alarm High (Hoog) wordt geactiveerd.



Afbeelding 3.41. Scherm High and Low alarms (Alarmen hoog en laag).

- Als u het alarm Low (Laag) wilt configureren, schuift u de knop On/Off (Aan/Off) naar de kant 'On' (Aan). Voer in het veld bij 'Low Alarm' (Laag alarm) de temperatuurwaarde in waarbij het alarm Low (Laag) wordt geactiveerd.
- 8. Nadat u de waarden voor de alarmen High (Hoog) en Low (Laag) hebt ingevoerd, raakt u 'Done' (Gereed) aan.
- 9. In het scherm Temperature Module Configuration (Module Temperatuur configureren) worden nu de ingevoerde waarden voor de alarmen High (Hoog) en Low (Laag) weergegeven. Als u de alarmen High High (Hoog-hoog)/Low Low (Laag-laag) wilt instellen voor temperatuur, raakt u het veld 'Low Low and High High alarms' (Alarmen Laag-laag en Hoog-hoog) aan.
- Het geopende scherm bevat aparte gedeeltes voor de alarmen High High (Hoog-hoog) en Low Low (Laag-laag). Schuif de knoppen On/Off (Aan/uit) in elk gedeelte naar de kant 'On' (Aan) om de alarmen in te schakelen.

11. Voor de alarmen High High (Hoog-hoog) en Low Low (Laaglaag) worden er velden met het label 'Interlock' en 'Hysteresis' (Hysterese) weergegeven. Raak 'Interlock' in een van de alarmvelden aan om de initiële voorwaarde in te stellen waarbij het alarm wordt geactiveerd, en 'Hysteresis' (Hysterese) om aan te geven welke actie na activering moet worden uitgevoerd.

\odot	Temperature High High and Low Low Alarms		
	High High Alarm	Off)	
	Interlock:	Temp1>0.00°C then Stop TCU	
	Hysteresis:	Resume if Temp1<0.00°C for 0s	
	Low Low Alarm	Off 🔵 🗕	
	Interlock:	Temp1<0.00°C then Stop TCU	
	Hysteresis:	Resume if Temp1>0.00°C for 0s	
		Cancel Done	

Afbeelding 3.42. Scherm High High and Low Low Alarms (Alarmen Hoog-hoog en Laag-laag).

- 12. In het geopende scherm (afbeelding 3.42) worden de gedeeltes 'Interlock' (Interlock) en 'Hysteresis' (Hysterese) getoond. In het gedeelte 'Interlock Equation' (Interlockvergelijking) voert u waarden in bij Control Set Point (Controle-instelpunt) en selecteert u een waarde bij Action Taken (Ondernomen actie) (Pause (Pauzeren) of Stop (Stoppen)).
- 13. Het gedeelte 'Hysteresis' (Hysterese) (onder het gedeelte 'Interlock') bepaalt wanneer processen worden hervat nadat het alarm is geactiveerd. Voer de tolerantie en de tijdsduur in gedurende welke het temperatuurniveau van het controleinstelpunt (plus de tolerantie) moet worden vastgehouden voordat wordt verdergegaan (in seconden). Wanneer u gereed bent, raakt u 'Done' (Gereed) aan.
- 14. Nadat u de interlock- en hystereseparameters voor de alarmen High High (Hoog-hoog) en Low Low (Laag-laag) hebt ingesteld, raakt u 'Done' (Gereed) aan. De Temperature Module Configuration (Module Temperatuur configureren) moet de waarden High High (Hoog-hoog) en/of Low Low (Laag-laag) tonen die u hebt opgegeven in het veld 'High High and Low Low Alarms' (Alarmen High High (Hoog-hoog) en Low Low (Laag-laag)).

- 15. Als u de temperatuur wilt kalibreren, raakt u het veld 'Calibration' (Kalibratie) aan. Volg de kalibratie-instructies uit hoofdstuk 4 — Kalibratieprocedures. U kunt de kalibratie ook overslaan en 'Done' (Gereed) aanraken om terug te gaan naar het startscherm.
- 16. Als u een TCU gebruikt, raadpleegt u de paragraaf over TCUconfiguratie en kalibratieparameters.

TCU-configuratie en kalibratieparameters

Het volgende schema illustreert de configuratie en kalibratieparameters voor TCU's uit de Lauda Integral T-serie, de Lauda Varicool-serie en de Neslab TF-serie. De blauwgroene vakken zijn items uit het menu Settings (Instellingen), de grijze vakken zijn parameters die verplicht moeten worden opgegeven.



Afbeelding 3.43. Configuratieparameters voor TCU uit de Neslab TF-serie.



Afbeelding 3.44. Configuratieparameters voor TCU uit de Lauda Varicool-serie.



Afbeelding 3.45. Configuratieparameters voor TCU uit de Lauda Integral T-serie.



Afbeelding 3.46. TCU-kalibratieprocedures voor Neslab TF-serie.



Afbeelding 3.47. TCU-kalibratieprocedures voor Lauda Varicool-serie.



Afbeelding 3.48. TCU-kalibratieprocedures voor Lauda Integral T-serie.

3.5.12 De module Liquid Pressure (Vloeistofdruk) configureren

Gebruik de volgende informatie en stappen om de module Liquid Pressure (Vloeistofdruk) te configureren. Elk van de onderstaande velden wordt geïllustreerd in afbeelding 3.49. **Opmerking:** U kunt 'Done' (Gereed) onder aan het scherm Configuration (Configuratie) te allen tijde aanraken tijdens de configuratie van de module. Uw invoer wordt opgeslagen en u kunt doorgaan waar u was gebleven wanneer u de moduleconfiguratie opnieuw opent.

Verplichte velden

• Units (Eenheden): Voer de eenheden in die u op de touchscreenconsole wilt gebruiken wanneer deze de vloeistofdruk weergeeft (mbar of psi).

Optionele velden

- Low and High alarms (Alarmen Laag en Hoog): Voer de vloeistofdrukwaarden in waarmee de vloeistofdrukalarmen High (Hoog) en Low (Laag) worden geactiveerd. Deze alarmen hebben geen invloed op lopende processen, maar waarschuwen gebruikers dat bepaalde omstandigheden buiten de gespecificeerde gebruiksparameters vallen.
- Low Low and High High alarms (Alarmen Laag-laag en Hoog-hoog): Vul de velden voor de interlockvergelijking en hysterese in om de voorwaarden in te stellen waarmee de vloeistofdrukalarmen High High (Hoog-hoog) en Low Low (Laaglaag) worden geactiveerd. Bij activering van deze alarmen klinkt een zoemer en worden lopende processen gepauzeerd of gestopt, afhankelijk van de gebruikersinstellingen. Voor de alarmen High High (Hoog-hoog) en Low Low (Laag-laag) moeten gebruikers waarden opgeven in de velden 'Interlock' en 'Hysteresis' (Hysterese). Deze waarden bepalen onder welke voorwaarden een alarm wordt geactiveerd, welke acties vervolgens moeten worden ondernomen en wanneer de functies worden hervat.
- Alarm designation (Alarmaanduiding): Nadat u de vloeistofdrukalarmen High High (Hoog-hoog) en/of Low Low (Laag-laag) hebt ingesteld, gebruikt u dit veld om te selecteren op welke pompen u deze alarmen wilt toepassen. Bijvoorbeeld voor Pomp 1 schakelt u het alarm High High (Hoog-hoog) in en het alarm Low Low (Laag-laag) uit.
- **Calibration (Kalibratie):** Configureer een nieuwe kalibratie voor vloeistofdruk met een éénpunts offset. Zie hoofdstuk 4 voor meer informatie over kalibratieprocedures.

 Zoek de module 'Liquid Pressure'(Vloeistofdruk) op het startscherm van de touchscreenconsole op en raak 'Configure' (Configureren) aan. Het scherm Liquid Pressure Module Configuration (Module Vloeistofdruk configureren) verschijnt (afbeelding 3.49). Raak het veld 'Units' (Eenheden) aan.

Liquid Pressure Module Configuration			
Liquid Pressure1			
Units:	PSI	•	
Low and High Alarms:	Lo Disabled-Hi Disab	•	
Low Low and High High Alarms:	LoLo Disabled-HiHi	~	
Alarm Designation (Pumps):	LoLo- None, HiHi- N	•	
Calibration:	Calibrate	•	
		Done	

Afbeelding 3.49. Scherm Liquid Pressure Module Configuration (Module Vloeistofdruk configureren).

- Selecteer mbar of psi in het geopende scherm. Raak 'Next' (Volgende) aan. Het scherm Liquid Pressure Module Configuration (Module Vloeistofdruk configureren) wordt opnieuw geopend. De geselecteerde eenheden worden weergegeven in het veld naast 'Units' (Eenheden).
- Als u de alarmen High (Hoog)/Low (Laag) wilt instellen voor vloeistofdruk, raakt u het veld 'Low and High alarms' (Alarmen Laag en Hoog) aan. Het geopende scherm bevat aparte gedeeltes voor de alarmen High (Hoog) en Low (Laag) (afbeelding 3.50). Als u het alarm High (Hoog) wilt configureren, schuift u de knop On/Off (Aan/Uit) naar de kant 'On' (Aan). Voer in het veld bij 'High Alarm' (Hoog alarm) de vloeistofdrukwaarde in waarbij het alarm High (Hoog) wordt geactiveerd. Raak 'Next' (Volgende) aan.



Afbeelding 3.50. Scherm Liquid Pressure High and Low Alarms (Alarmen Hoog en Laag voor vloeistofdruk).

- Als u het alarm Low (Laag) wilt configureren, schuift u de knop On/Off (Aan/Off) naar de kant 'On' (Aan). Voer in het veld bij 'Low Alarm' (Laag alarm) de vloeistofdrukwaarde in waarbij het alarm Low (Laag) wordt geactiveerd.
- 5. Nadat u de waarden voor de alarmen High (Hoog) en Low (Laag) hebt ingevoerd, raakt u 'Done' (Gereed) aan. In het scherm Liquid Pressure Module Configuration (Module Vloeistofdruk configureren) worden nu de ingevoerde waarden voor de alarmen High (Hoog) en Low (Laag) weergegeven.
- Als u de alarmen High High (Hoog-hoog)/Low Low (Laag-laag) wilt instellen voor vloeistofdruk, raakt u het veld 'Low Low and High High alarms' (Alarmen Laag-laag en Hoog-hoog) aan.
- 7. Voor de alarmen High High (Hoog-hoog) en Low Low (Laaglaag) worden er velden met het label 'Interlock' en 'Hysteresis' (Hysterese) weergegeven. Raak 'Interlock' in een van de alarmvelden aan om de initiële voorwaarde in te stellen waarbij het alarm wordt geactiveerd, of 'Hysteresis' (Hysterese) om aan te geven welke actie na activering moet worden uitgevoerd.
- In het geopende scherm worden de interlock- en hysteresegedeeltes getoond (afbeelding 3.51). In het gedeelte 'Agitation Interlock Equation' (Interlockvergelijking voor roeren) voert u waarden in bij Control Set Point (Controle-instelpunt) (in mbar of psi) en selecteert u een waarde bij Action Taken (Ondernomen actie) (Pause (Pauzeren) of Stop (Stoppen)).

$ \mathbf{\bullet} $	Liquid Pressure1 High High and Low Low Alarms		
	High High Alarm	Off 💿	
	Interlock:	Liquid Pressure1>0.00PSI then Stop	
	Hysteresis:	Resume if Liquid Pressure1<0.00PSI for 0s	
	Low Low Alarm	Off 🔵 🗕	
	Interlock:	Liquid Pressure1<0.00PSI then Stop	
	Hysteresis:	Resume if Liquid Pressure1>0.00PSI for 0s	
		Cancel Done	

Afbeelding 3.51. Scherm Liquid Pressure High High and Low Low Alarms (Alarmen Hoog-hoog en Laag-laag voor vloeistofdruk).

- Het gedeelte 'Hysteresis' (Hysterese) (onder het gedeelte Interlock) bepaalt de vloeistofdruk die moet worden bereikt (en hoe lang die vloeistofdruk moet worden aangehouden, in seconden) voordat wordt doorgegaan. Voer een waarde in bij Control Set Point (Controle-instelpunt) (in mbar of psi) en Seconds (Seconden). Wanneer u gereed bent, raakt u 'Done' (Gereed) aan.
- 10. Nadat u de interlock- en hystereseparameters voor de alarmen High High (Hoog-hoog) en Low Low (Laag-laag) hebt ingesteld, raakt u 'Done' (Gereed) aan. In het opnieuw geopende scherm Liquid Pressure Module Configuration (Module Vloeistofdruk configureren) worden nu de ingevoerde waarden voor de alarmen High High (Hoog-hoog) en Low Low (Laag-laag) weergegeven.
- 11. Als u een kalibratie voor vloeistofdruk wilt instellen, raakt u het veld 'Calibration' (Kalibratie) aan. Volg de kalibratie-instructies uit hoofdstuk 4 — Kalibratieprocedures. U kunt de kalibratie ook overslaan en 'Done' (Gereed) aanraken om terug te gaan naar het startscherm.

3.5.13 De module Auxiliary Output (Hulpuitgang) configureren

Gebruik de volgende informatie en stappen om de module Auxiliary Output (Hulpuitgang) te configureren. Elk van de onderstaande velden wordt geïllustreerd in afbeelding 3.52. **Opmerking:** U kunt 'Done' (Gereed) onder aan het scherm Configuration (Configuratie) te allen tijde aanraken tijdens de configuratie van de module. Uw invoer wordt opgeslagen en u kunt doorgaan waar u was gebleven wanneer u de moduleconfiguratie opnieuw opent.

Verplichte velden

- Sensor value to duplicate (Te dupliceren sensorwaarde): Selecteer het type sensorwaarde dat u wilt dupliceren. U kunt iedere beschikbare module op de touchscreenconsole selecteren, uitgezonderd Auxiliary Input (Hulpingang)
- Other required fields (Andere verplichte velden): Welke velden verplicht zijn voor de modules Auxiliary Output (Hulpuitgang) wisselt al naar gelang de module die u wilt dupliceren. Voordat u een module Auxiliary Output (Hulpuitgang) configureert, moet u eerst vertrouwd zijn met de module die u wilt dupliceren (paragrafen 3.5.2 tot 3.5.12). Deze paragrafen bevatten informatie over de verplichte en optionele velden in elke module.

Optionele velden

- Module name (Modulenaam): Voer een naam voor de module Auxiliary Ouput (Hulpuitgang) in. Het is raadzaam de naam in te voeren van de module die u dupliceert, samen met een nummer om de module Auxiliary Output (Hulpuitgang) aan te duiden. Als u bijvoorbeeld de module Conductivity (Geleidbaarheid) dupliceert, raden we aan de module Conductivity (Geleidbaarheid) de naam 'Geleidbaarheid 2' te geven om de modules van elkaar te onderscheiden.
- Zoek de module 'Auxiliary Output' (Hulpuitgang) op het startscherm van de touchscreenconsole op en raak 'Configure' (Configureren) aan. Het scherm Auxiliary Output Module Configuration (Module Hulpuitgang configureren) verschijnt (afbeelding 3.52). Raak het veld 'Sensor value to duplicate' (Te dupliceren sensorwaarde) aan.



Afbeelding 3.52. Scherm Auxiliary Output Module Configuration (Module Hulpuitgang configureren).

2. Raak in het geopende scherm (afbeelding 3.53) de module aan die u wilt dupliceren. Raak vervolgens 'Next' (Volgende) aan.

AuxOut4 Module Configuration		
	Which sensor would you like to duplicate to AuxOut4?	
	None	
	Agitation	
	BPC Pressure	
	Conductivity1	
	Conductivity2	
	Delta Pressure	
	Liquid Pressure1	
	Cancel Done	

Afbeelding 3.53. Module selecteren om te dupliceren.

- Het scherm Auxiliary Output Configuration (Hulpuitgang configureren) wordt opnieuw geopend met de verplichte en optionele velden van de module die u dupliceert. Zie de paragraaf over de configuratie van de module die u dupliceert (paragrafen 3.5.2 tot 3.5.12) voor informatie over alle bijbehorende velden.
- 4. Als u een naam wilt geven aan de module Auxiliary Output (Hulpuitgang), raak u het veld 'Module name' (Modulenaam) aan in het scherm Auxiliary Output Configuration (Hulpuitgang configureren). Geef de nieuwe modulenaam op met het geopende alfanumerieke toetsenbord. Raak 'Next' (Volgende) aan om terug te gaan naar het scherm Auxiliary Output Module Configuration (Module Hulpuitgang configureren).

3.5.14 De module Auxiliary Input (Hulpingang) configureren

Gebruik de volgende informatie en stappen om de module Auxiliary Input (Hulpingang) te configureren. Elk van de onderstaande velden wordt geïllustreerd in afbeelding 3.54. **Opmerking:** U kunt 'Done' (Gereed) onder aan het scherm Configuration (Configuratie) te allen tijde aanraken tijdens de configuratie van de module. Uw invoer wordt opgeslagen en u kunt doorgaan waar u was gebleven wanneer u de moduleconfiguratie opnieuw opent.

Verplichte velden

- Units (Eenheden): Voer de eenheden in die u op de touchscreenconsole wilt gebruiken voor de module Auxiliary Input (Hulpingang)
- Scaling (Schalen): Voer de schaling in die u op de touchscreenconsole wilt gebruiken voor het mA-signaal en de door u opgegeven eenheden.

Optionele velden

- Low and High alarms (Alarmen Laag en Hoog): Voer de waarden in waarmee de alarmen High (Hoog) en/of Low (Laag) worden geactiveerd voor de module Auxiliary Input (Hulpingang). Deze alarmen hebben geen invloed op lopende processen, maar waarschuwen gebruikers dat bepaalde omstandigheden buiten de gespecificeerde gebruiksparameters vallen.
- Low Low and High High alarms (Alarmen Laag-laag en Hoog-hoog): Vul de velden voor de interlockvergelijking en hysterese in om de voorwaarden in te stellen waarmee de hulpingangalarmen High High (Hoog-hoog) en/of Low Low (Laaglaag) worden geactiveerd. Bij activering van deze alarmen klinkt een zoemer en worden lopende processen gepauzeerd of gestopt, afhankelijk van de gebruikersinstellingen. Voor de alarmen High High (Hoog-hoog) en Low Low (Laag-laag) moeten gebruikers waarden opgeven in de velden 'Interlock' en 'Hysteresis' (Hysterese). Deze waarden bepalen onder welke voorwaarden een alarm wordt geactiveerd, welke acties vervolgens moeten worden ondernomen en wanneer de functies worden hervat.
- Alarm designation (Alarmaanduiding): Nadat u de alarmen High High (Hoog-hoog) en/of Low Low (Laag-laag) hebt ingesteld, gebruikt u dit veld om te selecteren op welke pompen u deze alarmen wilt toepassen. Bijvoorbeeld voor Pomp 1 schakelt u het alarm High High (Hoog-hoog) in en het alarm Low Low (Laag-laag) uit.
- Zoek de module 'Auxiliary Input' (Hulpingang) op het startscherm van de touchscreenconsole op en raak 'Configure' (Configureren) aan. Het scherm Auxiliary Input Module Configuration (Module Hulpingang configureren) verschijnt (zie afbeelding 3.54). Raak het veld 'Units' (Eenheden) aan.

AuxIn1 Module Configuration				
Units:	mA	•		
Scaling:	4.00mA - 4mA 20.00	•		
Low and high alarms:	Lo Disabled-Hi Disab	•		
Low low and high high alarms:	LoLo Disabled-HiHi	•		
Alarm designation (pumps):	LoLo- None, HiHi- N…	•		
		Done		

Afbeelding 3.54. Scherm Auxiliary Input Module Configuration (Module Hulpingang configureren).

- 2. Voer in het geopende scherm de eenheden in die u wilt gebruiken. Wanneer u gereed bent, raakt u 'Done' (Gereed) aan.
- 3. Het scherm Auxiliary Input Module Configuration (Module Hulpingang configureren) wordt opnieuw geopend. Raak het veld 'Scaling' (schaling) aan.
- 4. Voer in het geopende scherm de schaling in die u wilt gebruiken tussen het op de touchscreenconsole ontvangen mA-signaal en de eerder door u opgegeven eenheden. Wanneer u gereed bent, raakt u 'Done' (Gereed) aan.
- Het scherm Auxiliary Input Module Configuration (Module Hulpingang configureren) wordt opnieuw geopend. Als u de alarmen High (Hoog)/Low (Laag) wilt instellen voor de module Auxiliary Input (Hulpingang), raakt u het veld 'Low en High alarms' (Alarmen Laag en Hoog) aan.
- Het geopende scherm bevat aparte gedeeltes voor de alarmen High (Hoog) en Low (Laag). Als u het alarm High (Hoog) wilt configureren, schuift u de knop On/Off (Aan/Uit) naar de kant 'On' (Aan). Voer in het veld bij 'High Alarm' (Hoog alarm) de waarde in waarbij het alarm High (Hoog) wordt geactiveerd. Raak 'Next' (Volgende) aan.
- Als u het alarm Low (Laag) wilt configureren, schuift u de knop On/Off (Aan/Off) naar de kant 'On' (Aan). Voer in het veld bij 'Low Alarm' (Laag alarm) de waarde in waarbij het alarm Low (Laag) wordt geactiveerd.

- Nadat u de waarden voor de alarmen High (Hoog) en Low (Laag) hebt ingevoerd, raakt u 'Done' (Gereed) aan. In het scherm Auxiliary Input Module Configuration (Module Hulpingang configureren) worden nu de ingevoerde waarden voor de alarmen High (Hoog) en Low (Laag) weergegeven.
- Als u de alarmen High High (Hoog-hoog)/Low Low (Laag-laag) wilt instellen voor de module Auxiliary Input (Hulpingang), raakt u het veld 'Low low en High high alarms' (Alarmen Laag-laag en Hooghoog) aan.
- 10. Voor de alarmen High High (Hoog-hoog) en Low Low (Laaglaag) worden er velden met het label 'Interlock' en 'Hysteresis' (Hysterese) weergegeven. Raak 'Interlock' in een van de alarmvelden aan om de initiële voorwaarde in te stellen waarbij het alarm wordt geactiveerd, of 'Hysteresis' (Hysterese) om aan te geven welke actie na activering moet worden uitgevoerd.
- 11. In het geopende scherm worden de interlock- en hysteresegedeeltes getoond. In het gedeelte 'Interlock Equation' (Interlockvergelijking) voert u waarden in bij Control Set Point (Controle-instelpunt) en selecteert u een waarde bij Action Taken (Ondernomen actie) (Pause (Pauzeren) of Stop (Stoppen)).
- 12. Het gedeelte 'Hysteresis' (Hysterese) (onder het gedeelte Interlock) bepaalt de waarde die moet worden bereikt (en hoe lang die waarde moet worden aangehouden, in seconden) voordat wordt doorgegaan. Voer een waarde in bij Control Set Point (Controleinstelpunt) en Seconds (Seconden). Wanneer u gereed bent, raakt u 'Done' (Gereed) aan.
- 13. Nadat u de interlock- en hystereseparameters voor de alarmen High High (Hoog-hoog) en Low Low (Laag-laag) hebt ingesteld, raakt u 'Done' (Gereed) aan. In het opnieuw geopende scherm Auxiliary Input Module Configuration (Module Hulpingang configureren) worden nu de ingevoerde waarden voor de alarmen High High (Hoog-hoog) en/of Low Low (Laag-laag) weergegeven.

4

Kalibratieprocedures

Inhoudsopgave

4.1

- Pompen kalibreren
- 4.2 pH kalibreren
- 4.3 Geleidbaarheid kalibreren
- 4.4 Roersnelheid kalibreren
- 4.5 Massa kalibreren
- 4.6 BPC-luchtdruk kalibreren
- 4.7 Vloeistofdruk kalibreren
- 4.8 Temperatuur kalibreren

4.1 Pompen kalibreren

- 1. Als u nog geen actieve pompkalibratie hebt ingesteld, wordt u nu in een scherm gevraagd een nieuwe kalibratie te maken. Raak de knop 'Calibrate now' (Nu kalibreren) aan.
- Vul de velden in voor het slangtype, de binnendiameter van slangen, het vloeistoftype, de eenheden en de vloeistofdichtheid. U kunt desgewenst de pompkalibratie ook een naam geven en opmerkingen toevoegen. Raak 'Next' (Volgende) aan.
- 3. Gebruikers kunnen kiezen of ze massa of volume willen gebruiken om pompen te kalibreren. Voor massakalibratie hebt u een beker en een schaal nodig. Een cilinder met maataanduiding is vereist voor volumekalibratie. Voor de optie Volume moeten gebruikers de eenheden opgeven (ml of I), voor de optie Mass (Massa) alleen grammen. Voor beide opties moeten gebruikers RPM-limieten (maximale en minimale snelheden) en vloeistofdichtheid opgeven.
- 4. Nadat u de verplichte velden hebt ingevuld, kunt u 'Prime Set Up' (Primen configureren) of 'Skip Prime' (Primen overslaan) opgeven. Als u 'Prime Set Up' (Primen configureren) selecteert, wordt u in een scherm gevraagd de beker of een cilinder met maataanduiding op een schaal te configureren.
 - a. Nadat u 'Next' (Volgende) hebt aangeraakt, wordt u in een tweede scherm gevraagd slangen in de pomp te laden en de slangen vervolgens in de beker of cilinder met maataanduiding te plaatsen.
 - B. Raak 'Next' (Volgende) nogmaals aan. Houd de knop 'Prime' (Primen) ingedrukt totdat het instelpunt voor primen op het scherm wordt weergegeven.
 - c. Selecteer 'Next' (Volgende). U wordt doorgestuurd naar het scherm 'Calibration Total Time Input' (Totale kalibratietijd invoeren) en wordt gevraagd de totale tijd voor pompkalibratie in te voeren (in seconden).

Opmerking: Als u 'Skip Prime' (Primen overslaan) kiest, wordt u rechtstreeks doorgestuurd naar het scherm 'Calibration Total Time Input' (Totale kalibratietijd invoeren). Voer de totale tijd voor pompkalibratie in seconden in.

5. Er zijn vier kalibratiepunten voor de volume- en massaopties, en vier stappen voor elk kalibratiepunt. Herhaal de volgende stappen voor elk van de vier kalibratiepunten:

- a. Kies om de beker/cilinder met maataanduiding te legen en de schaal na elk kalibratiepunt op nul te stellen, of om de beker/ cilinder met maataanduiding te vullen en de schaal na elk kalibratiepunt niet op nul te stellen.
- b. Configureer een lege beker/cilinder met maataanduiding en een schaal. Tarreer de weegschaal, zodat alleen het gewicht van de vloeistof wordt gemeten en niet het gewicht van de vloeistof en de beker. Raak 'Next' (Volgende) aan.
- c. Raak de knop 'Deliver Bolus' (Bolus toedienen) aan om een bolus toe te voegen aan de beker/cilinder met maataanduiding. Voor elk kalibratiepunt voert de pomp een ander percentage van het beschikbare RPM-uitvoerbereik uit (van laag naar hoog) gedurende een verschillende tijdsduur.
- d. Nadat de bolus is toegediend, verschijnt de melding 'Delivery Complete' (Toediening voltooid). Raak 'Next' (Volgende) aan. Voer de totale hoeveelheid vloeistof in die is gepompt tijdens het kalibratiepunt (1, 2, 3 of 4) en raak 'Calibrate' (Kalibreren) aan. Het scherm voor het volgende kalibratiepunt wordt geopend.
- Nadat alle vier de kalibratiepunten zijn voltooid, raakt u 'Done' (Gereed) rechtsonder op het scherm aan om terug te gaan naar het scherm Pump Settings (Pompinstellingen).

4.2 pH kalibreren

- Als u nog geen actieve pH-kalibratie hebt ingesteld, wordt u nu in een scherm gevraagd een nieuwe kalibratie te maken. Raak de knop 'Create New Calibration' (Nieuwe kalibratie maken) aan. Als u al een actieve pH-kalibratie hebt gemaakt, raakt u pH-module op het startscherm van de touchscreenconsole aan om het scherm 'pH Settings' (pH-instellingen) te openen. Raak de knop 'Calibration' (Kalibreren) aan.
- 2. Selecteer eerst welke pH-sensor u kalibreren. Kies vervolgens het type kalibratie dat u wilt gebruiken (geef helling en offset, offsetkalibratie, 2-punts of 3-punts op).
- 3. Volg de onderstaande stappen om de helling en offset te kalibreren:
 - Nadat u de knop 'Enter slope and offset' (Helling en offset opgeven) hebt aangeraakt, wordt de gebruiker in een scherm gevraagd de helling (pH/mV) en offset (mV) op te geven.
 - b. Raak rechtsonder in het scherm 'Done' (Gereed) aan om terug te gaan naar het scherm pH Settings (pH-instellingen).

- 4. Volg de onderstaande stappen om de offset te kalibreren:
 - a. Nadat u de knop 'Offset Calibration' (Offsetkalibratie) hebt aangeraakt, wordt er een scherm geopend met de stabiliserende pH-sensorwaarde voor kalibratiepunt 1. Nadat de waarde is gestabiliseerd, wordt het veld 'Buffer Solution Value' (Waarde bufferoplossing) automatisch ingevuld en precies afgestemd op de bufferwaarde. U kunt de waarde desgewenst ook wijzigen. Raak 'Calibrate' (Kalibreren) aan.
 - b. Raak rechtsonder in het scherm 'Done' (Gereed) aan om terug te gaan naar het scherm pH Settings (pH-instellingen).
- Volg de onderstaande stappen voor 2-punts kalibratie.
 Opmerking: Voor driepunts kalibratie worden dezelfde stappen gevolgd, maar worden stappen 'd' en 'e' voor het derde kalibratiepunt herhaald.
 - a. Nadat u de knop '2-Point Calibration' (2-punts kalibratie) hebt aangeraakt, verschijnt er een scherm met het verzoek om een temperatuurcompensatie voor de kalibratie te selecteren. De verwachte buffertemperatuur wordt gebruikt als er geen RTD aanwezig is. Indien aanwezig kunt u de RTD ook gebruiken door de schuifknop in te stellen om deze in te schakelen. Raak 'Next' (Volgende) aan.
 - b. In het geopende scherm wordt de gebruiker gevraagd de pH-sensor (en RTD, indien geselecteerd) te plaatsen in de eerste bufferoplossing voor kalibratiepunt 1. Raak 'Next' (Volgende) aan.
 - c. Het volgende scherm toont de stabiliserende pHsensorwaarde voor kalibratiepunt 1. Nadat de waarde is gestabiliseerd, wordt het veld 'pH Value' (pH-waarde) automatisch ingevuld en precies afgestemd op de bufferwaarde. U kunt de waarde desgewenst ook wijzigen. Raak 'Calibrate' (Kalibreren) aan.
 - In het geopende scherm wordt de gebruiker gevraagd de pH-sensor (en RTD, indien geselecteerd) te plaatsen in de tweede bufferoplossing voor kalibratiepunt 2. Raak 'Next' (Volgende) aan.
 - e. Het volgende scherm toont de stabiliserende pHsensorwaarde voor kalibratiepunt 2. Nadat de waarde is gestabiliseerd, wordt het veld 'pH Value' (pH-waarde) automatisch ingevuld en precies afgestemd op de bufferwaarde. U kunt de waarde desgewenst ook wijzigen. Raak 'Calibrate' (Kalibreren) aan.
 - f. Het volgende scherm meldt dat de pH-kalibratie gereed is. De gebruiker wordt gevraagd de gewenste temperatuurcompensatie in het volgende scherm bij te werken. Raak 'Next' (Volgende) aan.

- g. Stel de temperatuurcompensatie in die moet worden gebruikt als de sensor is ingesteld als actief. De verwachte buffertemperatuur wordt gebruikt als er geen RTD aanwezig is. Indien aanwezig kunt u de RTD ook gebruiken door de schuifknop in te stellen om deze in te schakelen. Raak 'Done' (Gereed) aan.
- h. Het scherm 'pH Calibration Complete' (pH-kalibratie voltooid) verschijnt. Raak rechtsonder in het scherm 'Done' (Gereed) aan om terug te gaan naar het scherm pH Settings (pH-instellingen).

4.3 Geleidbaarheid kalibreren

- Als u nog geen actieve geleidbaarheidskalibratie hebt ingesteld, wordt u in een scherm gevraagd een nieuwe kalibratie te maken. Raak de knop 'Create New Calibration' (Nieuwe kalibratie maken) aan. Als u al een actieve geleidbaarheidskalibratie hebt gemaakt, raakt u module Conductivity (Geleidbaarheid) op het startscherm van de touchscreenconsole aan om het scherm 'Conductivity Settings' (Geleidbaarheidsinstellingen) te openen. Raak de knop 'Calibration' (Kalibreren) aan.
- In het geopende scherm wordt de gebruiker gevraagd een geleidbaarheidssensor voor kalibratie te selecteren (Conductivity1 of Conductivity2). Nadat u een geleidbaarheidssensor hebt geselecteerd, wordt u gevraagd welke kalibratiemethode u wilt gebruiken (1-punts, 2-punts of 3-punts kalibratie).
- 3. Volg de onderstaande stappen voor 1-punts kalibratie:
 - a. Voer in het geopende scherm de celconstante van de geleidbaarheidssensor in. Raak vervolgens 'Next' (Volgende) aan.
 - b. In het geopende scherm wordt de gebruiker gevraagd de geleidbaarheidssensor in een bufferoplossing te plaatsen.
 Raak 'Next' (Volgende) aan.
 - c. Het volgende scherm toont de stabiliserende geleidbaarheidssensorwaarde voor kalibratiepunt 1 van 1 (in µS/cm). Nadat de waarde is gestabiliseerd, wordt het veld 'Conductivity Value' (Geleidbaarheidswaarde) automatisch ingevuld en precies afgestemd op de bufferwaarde. U kunt de waarde desgewenst ook wijzigen. Raak 'Calibrate' (Kalibreren) aan.

- d. Wanneer het scherm 'Conductivity Calibration Complete' (Geleidbaarheidskalibratie voltooid) verschijnt, is de kalibratie voltooid. Raak rechtsonder in het scherm 'Done' (Gereed) aan om terug te gaan naar het scherm Conductivity Settings (Geleidbaarheidsinstellingen).
- Volg de onderstaande stappen voor 2-punts kalibratie.
 Opmerking: Voor 3-punts kalibratie worden dezelfde stappen gevolgd, maar worden stappen 'd' en 'e' voor het derde kalibratiepunt herhaald.
 - a. Voer in het geopende scherm de celconstante van de geleidbaarheidssensor in. Raak vervolgens 'Next' (Volgende) aan.
 - b. In het geopende scherm wordt de gebruiker gevraagd de geleidbaarheidssensor in de eerste bufferoplossing te plaatsen. Raak 'Next' (Volgende) aan.
 - c. Het volgende scherm toont de stabiliserende geleidbaarheidssensorwaarde voor kalibratiepunt 1 (in μS/cm). Nadat de waarde is gestabiliseerd, wordt het veld 'Conductivity Value' (Geleidbaarheidswaarde) automatisch ingevuld en precies afgestemd op de bufferwaarde. U kunt de waarde desgewenst ook wijzigen. Raak 'Calibrate' (Kalibreren) aan.
 - In het geopende scherm wordt de gebruiker gevraagd de geleidbaarheidssensor in de volgende bufferoplossing te plaatsen. Raak 'Next' (Volgende) aan.
 - e. Het volgende scherm toont de stabiliserende geleidbaarheidssensorwaarde voor het kalibratiepunt (in μS/cm). Nadat de waarde is gestabiliseerd, wordt het veld 'Conductivity Value' (Geleidbaarheidswaarde) automatisch ingevuld en precies afgestemd op de bufferwaarde. U kunt de waarde desgewenst ook wijzigen. Raak 'Calibrate' (Kalibreren) aan.
 - f. Wanneer het scherm 'Conductivity Calibration Complete' (Geleidbaarheidskalibratie voltooid) verschijnt, is de kalibratie voltooid. Raak rechtsonder in het scherm 'Done' (Gereed) aan om terug te gaan naar het scherm Conductivity Settings (Geleidbaarheidsinstellingen).

4.4 Roersnelheid kalibreren

 Raak de module Agitation (Roeren) in het startscherm van de touchscreenconsole aan om het scherm 'Agitation Settings' (Roer-instellingen) te openen. Raak de knop 'Calibration' (Kalibreren) aan.

- Er verschijnt een scherm waarin de gebruiker wordt gevraagd de werkelijke motorsnelheid (in RPM) te meten met een tachometer en de snelheid in het betreffende veld in te voeren. Nadat u de werkelijke motorsnelheid hebt ingevoerd, raakt u 'Done' (Gereed) aan.
- Het scherm '1-point offset adjustment complete' (Afstelling 1-punts kalibratie voltooid) verschijnt wanneer de kalibratie voltooid is. Raak rechtsonder in het scherm 'Done' (Gereed) aan om terug te gaan naar het scherm 'Agitation Settings' (Roer-instellingen).

4.5 Massa kalibreren

- Raak de module Mass (Massa) in het startscherm van de touchscreenconsole aan om het scherm 'Mass Settings' (Massa-instellingen) te openen. Raak de knop 'Calibration' (Kalibreren) aan.
- Selecteer het type kalibratie dat u wilt gebruiken (1-punts, 2-punts of 3-punts). **Opmerking:** Hierna vindt u de stappen van het proces voor 2-punts kalibratie. Voor 1-punts en 3-punts kalibratie worden dezelfde stappen gevolgd, met een paar kleine verschillen: 1-punts kalibratie heeft geen stap 'b', terwijl voor 3-punts kalibratie stap 'b' wordt herhaald voor het derde kalibratiepunt.
- 3. Volg de onderstaande stappen voor 2-punts kalibratie:
 - a. Nadat u de knop '2-point calibration' (2-punts kalibratie) hebt aangeraakt, verschijnt er voor 'Calibration Point 1 of 2' (Knoppen 1 van 2) een scherm waarin de gebruiker wordt gevraagd de eerste massasensorwaarde op te geven in kg. Plaats een gewicht in het vat, voer de sensorwaarde op het scherm in en raak 'Calibrate' (Kalibreren) aan.
 - b. Er verschijnt een scherm voor 'Calibration Point 2 of 2' (Kalibratiepunt 2 van 2). Plaats een tweede gewicht dat zwaarder is dan dat uit de eerste stap, in het vat. Voer de sensorwaarde van de totale massa in kg in en raak vervolgens 'Calibrate' (Kalibreren) aan.
 - c. Wanneer het scherm 'Mass Calibration Complete' (Geleidbaarheidskalibratie voltooid) verschijnt, is de kalibratie voltooid. Raak rechtsonder in het scherm 'Done' (Gereed) aan om terug te gaan naar het scherm 'Mass Settings' (Massa-instellingen).

4.6 BPC-luchtdruk kalibreren

- Als u nog geen actieve BPC-drukkalibratie hebt ingesteld, wordt u in een scherm gevraagd een nieuwe kalibratie te maken. Raak de knop 'Create New Calibration' (Nieuwe kalibratie maken) aan. Als u al een actieve BPC-drukkalibratie hebt gemaakt, raakt u module BPC Pressure (BPC-druk) op het startscherm van de touchscreenconsole aan om het scherm 'BPC Pressure Settings' (BPC-drukinstellingen) te openen. Raak de knop 'Calibration' (Kalibreren) aan.
- 2. In een scherm wordt de gebruiker gevraagd een kalibratiemethode te kiezen: nulstellen/tarreren, een 1-punts offsetwaarde invoeren of offset wissen. Selecteer 'Enter a 1-point offset value' (1-punts offsetwaarde invoeren').
- 3. In het volgende scherm wordt de gebruiker gevraagd de BPCsensorwaarde in psi op te geven. Raak 'Calibrate' (Kalibreren) aan nadat u de BPC-druk in psi hebt opgegeven.
- 4. Wanneer het scherm 'BPC Pressure Calibration Complete' (BPCdrukkalibratie voltooid) verschijnt, is de kalibratie voltooid. Raak rechtsonder in het scherm 'Done' (Gereed) aan om terug te gaan naar het scherm 'BPC Pressure Settings' (BPC-drukinstellingen).

4.7 Vloeistofdruk kalibreren

- Als u nog geen actieve vloeistofdrukkalibratie hebt ingesteld, wordt u in een scherm gevraagd een nieuwe kalibratie te maken. Raak de knop 'Create New Calibration' (Nieuwe kalibratie maken) aan. Als u al een actieve vloeistofdrukkalibratie hebt gemaakt, raakt u module Liquid Pressure (Vloeistofdruk) op het startscherm van de touchscreenconsole aan om het scherm Liquid Pressure Settings' (vloeistofdrukinstellingen) te openen. Raak de knop 'Calibration' (Kalibreren) aan.
- In een scherm wordt de gebruiker gevraagd een kalibratiemethode te kiezen: vloeistofdruk1 nulstellen/tarreren, vloeistofdruk2 nulstellen/tarreren, een 1-punts offsetwaarde invoeren of offset wissen. Selecteer 'Enter a 1-point offset value' (1-punts offsetwaarde invoeren').
- 3. Selecteer in het volgende scherm een vloeistofdruksensor die u voor kalibratie wilt gebruiken.

- In het volgende scherm wordt de gebruiker gevraagd de sensorwaarde voor vloeistofdruk in psi op te geven. Raak 'Calibrate' (Kalibreren) aan nadat u de vloeistofdruk in psi hebt opgegeven.
- Wanneer het scherm 'Liquid Pressure Calibration Complete' (Vloeistofdrukkalibratie voltooid) verschijnt, is de kalibratie voltooid. Raak rechtsonder in het scherm 'Done' (Gereed) aan om terug te gaan naar het scherm 'Liquid Pressure Settings' (Vloeistofdrukinstellingen).

4.8 Temperatuur kalibreren

- Als u nog geen actieve temperatuurkalibratie hebt ingesteld, wordt u in een scherm gevraagd een nieuwe kalibratie te maken. Raak de knop 'Create New Calibration' (Nieuwe kalibratie maken) aan. Als u al een actieve temperatuurkalibratie hebt gemaakt, raakt u module Temperature (Temperatuur) op het startscherm van de touchscreenconsole aan om het scherm 'Temperature Settings' (Temperatuurinstellingen) te openen. Raak de knop 'Calibration' (Kalibreren) aan.
- 2. In het geopende scherm wordt de gebruiker gevraagd een temperatuursensor voor kalibratie te selecteren.
- In het volgende scherm wordt de gebruiker gevraagd een kalibratiemethode te selecteren. Selecteer '1-Point Calibration' (1-punts kalibratie). **Opmerking:** De hier gegeven instructies illustreren 1-punts kalibratie. Herhaal de onderstaande stappen voor 2-punts en 3-punts kalibratie.
- In het volgende scherm wordt de gebruiker gevraagd de temperatuursensorwaarde op te geven (in°C). Raak 'Calibrate' (Kalibreren) aan nadat u de temperatuursensorwaarde hebt opgegeven.
- 5. Wanneer het scherm 'Temperature Calibration Complete' (Temperatuurkalibratie voltooid) verschijnt, is de kalibratie voltooid. Raak rechtsonder in het scherm 'Done' (Gereed) aan om terug te gaan naar het scherm 'Temperature Settings' (Temperatuurinstellingen).

6

BPC laden en sonde inbrengen

Inhoudsopgave

- 5.1 Richtlijnen voor algemeen gebruik
- 5.2 De BPC laden
- 5.3 De recirculatielijn configureren
- 5.4 De BPC met lucht vullen
- 5.5 De aandrijfas inbrengen
- 5.6 BPC's laden voor open-top mengen
- 5.7 Sondeaansluitingen maken
- 5.8 Sensoren en sondes inbrengen
- 5.9 De S.U.M. tarreren

5.1 Richtlijnen voor algemeen gebruik

5.1.1 BPC voorbereiden en configureren

Maak uzelf vertrouwd met de S.U.M. BPC en hardware voordat de BPC wordt geladen. De voorzijde van de BPC is ter referentie aangewezen als het paneel met de lagerpoort. De lagerpoort moet wijzen naar de ontvanger van de lagerpoort. Deze bevindt zich onder de motorsteun wanneer de BPC wordt geladen. Controleer bovendien of het volume van de BPC die wordt geladen, overeenkomt met het volume van de S.U.M.

5.1.2 Hanteringsinstructies BPC

Gebruik geen schaar of andere scherpe voorwerpen om de buitenste plastic zakken van de BPC's te openen. Wanneer u een BPC in de buitenste ondersteuningscontainer plaatst, sleep de BPC dan niet over hoeken of scherpe voorwerpen. Til de BPC niet aan de hoeken of bovenste naden. Rol de slangen bovenop de BPC voorzichtig op voor opslag om te voorkomen dat ze doorboord worden door kabelbinders of kabelklemmen. Gebruik beschermingsmateriaal tussen de slangen en de BPC voor opslag en transport.

5.1.3 Werkvolume

ledere S.U.M. is ontworpen voor een specifiek werkvolumebereik. Het minimale werkvolume en het nominale werkvolume worden vermeld in hoofdstuk 8—Specificaties en informatie over onderdelen. Voor normaal gebruik mogen de werkelijke werkvolumes de aangegeven nominale werkvolumes niet overschrijden.

5.1.4 Vloeistofoverdracht

De S.U.M. BPC is ontworpen met thermoplastische slangen, snelaansluitingen en drievoudige klemmen om vloeistof toe te voegen aan en te onttrekken uit de BPC. Voor toepassingen met vloeistof op vloeistof kan een steriele omgeving worden behouden zolang alle verbindingen op een aseptische manier worden gemaakt. Om de steriliteit van het mengsysteem te behouden, moeten extra leidingen (snelaansluitingen en drieklem) worden aangesloten onder een laminaire luchtstroomkap of met een steriele buizenlasser.
5.1.5 Toediening

Het roermechanisme mag niet worden bediend wanneer de volumes kleiner zijn dan het vermelde minimumvolume. De S.U.M. heeft een afvoerlijn waarmee vloeistof verwijderd kan worden met een peristaltische pomp. De afvoer bevindt zich aan de onderkant van de S.U.M. BPC, zodat er vrijwel geen verstoppingen plaatsvinden. De onderste afvoerlijn kan aangesloten worden met behulp van de meegeleverde snelaansluiting van 12,7 mm (0,5 inch). Schakel het roeren tijdens het aftappen uit wanneer het minimale werkvolume bijna bereikt is.

5.2 De BPC laden

Volg de onderstaande stappen om een BPC in de buitenste ondersteuningscontainer te laden.

Opmerking: Het is van cruciaal belang dat deze procedures worden gevolgd om de S.U.M. succesvol te kunnen gebruiken. Voor grotere systemen (500 tot 2.000 l) raden we aan dat twee gebruikers de BPC in de buitenste ondersteuningscontainer plaatsen. Voor grotere systemen is tevens toegang tot een ladder of een ander verhoogd platform vereist.

1. Als u een S.U.M. van 500 l, 1.000 l of 2.000 l gebruikt, opent u eerst de deur van de buitenste ondersteuningscontainer (afbeelding 5.1).



Afbeelding 5.1. De deur op een S.U.M. van 500 l openen.

2. Haal de BPC uit de dubbele plastic beschermingszakken (afbeelding 5.2). Let op dat u niet in de BPC knipt of snijdt wanneer u de plastic zak opent.



Afbeelding 5.2. De plastic zak openen.

- Nadat u de BPC uit de dubbele plastic zak hebt verwijderd, inspecteert u de BPC visueel op beschadiging. Als de BPC is beschadigd, neemt u onmiddellijk contact op met uw verkoopvertegenwoordiger.
- 4. Sluit alle klemmen van de BPC-leiding (afbeelding 5.3) Zorg dat de klem van de afvoerleiding zich zo dicht mogelijk bij de BPC-poort bevindt en volledig is gesloten.



Afbeelding 5.3. De klem van de leidingset sluiten.

 Richt de BPC met de lagerpoort omhoog naar de motoraandrijving en met de sondepoorten (indien aanwezig) naar de onderkant van de onderste toegangsuitsparing. Laad bij S.U.M.'s van 500 l, 1.000 l of 2.000 l de BPC door de open deur van de buitenste ondersteuningscontainer. Laad bij S.U.M.'s van 50 l, 100 l of 200 l de BPC via de bovenkant van de buitenste ondersteuningscontainer. Voer de afvoerleiding door de onderste bodemplaat (afbeelding 5.4). De bodemplaat kan worden verwijderd om het laden van deze onderste leidingset te vergemakkelijken. Positioneer de BPC vervolgens in de tank. Bij grotere units sluit u de deur op de buitenste ondersteuningscontainer.



Afbeelding 5.4. De afvoerleiding door de tank voeren.

7. Als u een BPC met sondepoorten gebruikt, moet u controleren of alle poortklemmen op de aansluitleidingen van de poort zijn gesloten en zich zo dicht mogelijk bij de behuizing van de BPC bevinden (afbeelding 5.5).



Afbeelding 5.5. De locatie van de poortklemmen controleren.

8. Open de vergrendeling van de lagerpoort bij de voorkant van het motorblok. Controleer of de zwarte bumpers aanwezig zijn op de achterkant van het motorsteunblok en aan de binnenkant van de vergrendeling (afbeelding 5.6). Breng de lagerpoort aan in de ontvanger. Sluit vervolgens de vergrendeling en klem deze vast (afbeelding 5.7).



Afbeelding 5.6. De lagerpoort in de ontvanger inbrengen.



Afbeelding 5.7. De ontvanger van de lagerpot sluiten en vergrendelen.

9. Gebruik de vier onderste uitsparingen in de basis van de buitenste ondersteuningscontainer om de positioneringstabs op de BPC uit te lijnen op de hardware (afbeelding 5.8).



Afbeelding 5.8. De BPC-positioneringstabs uitlijnen met pennen.

- 10. Bevestig de lipjes op de BPC aan elk van de vier onderste hoekpennen op de buitenste ondersteuningscontainer. Uitsparingen in de achterkant van de tank vergemakkelijken de bevestiging van de lipjes. Bij units van 500 l en 1.000 l kunt u de bevestiging van de lipjes vergemakkelijken door de beugel omhoog te trekken over de uitsparing in de achterkant van de tank en deze te verwijderen. Bij units van 2.000 l is een deur in de achterkant van de tank voorzien om de bevestiging van de positioneringslipjes aan de pennen te vergemakkelijken.
- 11. Wanneer de BPC volledig is aangebracht, moet u zorgen dat de afvoerleiding volledig door de onderste poort wordt geleid en dat de lagerpoort op de BPC zich in de juiste positie bevindt (wijzend naar de ontvangerpoort).
- 12. Zodra de lagerpoort volledig vastzit, moet u controleren of de film niet te strak rond de lagerpoort wordt getrokken. Het betreffende gebied bevindt zich op het voorste BPC-paneel, onder de lagerpoort (afbeeldingen 5.9 en 5.10). Om juiste inbreng van de aandrijfas te waarborgen en de spanning op de lagerpoort te verminderen, trekt u het voorpaneel van de BPC iets omhoog om extra film aan het betreffende gebied toe te voegen.



Afbeelding 5.9. Juist geladen BPC



Afbeelding 5.10. Onjuist geladen BPC

- 13. Als de gebruikte BPC sondepoorten heeft, lijnt u de rij sondepoorten uit via het toegangsvenster.
- 14. Als u systeem knijpventielen heeft, leidt u de slangen van de BPC door de knijpventielen. Knijpventielen kunnen worden opgeborgen op een optionele hanger aangebracht op de buitenste ondersteuningscontainer.

5.3 De recirculatielijn configureren

U kunt desgewenst een optionele recirculatie- of monsternamelus configureren Volg de onderstaande stappen om de recirculatielijn aan te sluiten.

1. Voer de afvoerleiding door een peristaltische pomp (afbeelding 5.11).



Afbeelding 5.11. Afvoerleiding in peristaltische pomp.

 Voeg bij niet-steriele toepassingen de snelaansluitingen van de recirculatie- of monsternamelijn en de afvoerleiding samen om de recirculatielus te vormen. Om deze verbinding te maken, opent u eerst de tamper-proof plastic zak die de snelaansluitingen op de recirculatie-/monsternamelijn en de afvoerleidingen bedekt (afvoerleiding 5.12).



Afbeelding 5.12. Plastic zak verwijderen.

 Haal de doppen van de snelaansluitingen. Lijn de snelaansluitingen van de recirculatie- en afvoerlijn uit en verbind ze (afbeelding 5.13).
 Opmerking: Bij steriele toepassingen moet deze verbinding worden aangebracht onder een laminaire luchtstroomkap of met gebruik van een steriele slanglasser.



Afbeelding 5.13. Leiding aansluiten.

4. Zorg dat de klemmen op de recirculatielijn zijn gesloten voordat u de BPC vult.

5. Wanneer u klaar bent om de recirculatielus te gebruiken, opent u alle klemmen op beide leidingensets. Afbeelding 5.14 toont een voltooide recirculatielus.



Afbeelding 5.14. Volledige recirculatielus.

5.4 De BPC met lucht vullen

Voordat u begint de BPC met lucht te vullen, moet u controleren of de BPC-drukmodule is geconfigureerd op de touchscreenconsole en of het S.U.M.-systeem is aangesloten op de luchttoevoer van uw faciliteit.

- 1. Raak de module 'BPC Pressure' (BPC-druk) op het startscherm van de touchscreenconsole aan. Selecteer 'Calibration' (Kalibratie)
- 2. Raak 'Zero/Tare BPC Pressure Sensor' (BPC-druksensor op nul stellen/tarreren) aan en voer 0 (nul) psi in het veld in
- 3. Raak 'Close' (Sluiten) onder aan het scherm BPC Pressure Module (BPC-drukmodule) aan om terug te gaan naar het startscherm.
- 4. Raak 'Start' (Starten) aan om te beginnen met het vullen met lucht. **Opmerking:** De touchscreenconsole heeft een ingeprogrammeerd instelpunt van 0,06 psi.

5.5 De aandrijfas inbrengen

Voordat u de aandrijfas inbrengt, moet u zich vertrouwd maken met de onderstaande illustratie van de mengeenheid (afbeelding 5.15). **Opmerking:** Wanneer u de aandrijfas inbrengt of verwijdert, moet u voorkomen dat de as tegen het plafond of plafondlampen stoot. Zie hoofdstuk 8 voor de vereiste plafondhoogte om de aandrijfas aan te brengen.



Afbeelding 5.15. Mengeenheid

- 1. Voordat u de aandrijfas laadt, moet u controleren of de BPC is gevuld met lucht (paragraaf 5.4).
- Verwijder de veiligheidsdeksel van de motor (afbeelding 5.16). Gebruik de twee bijgeleverde sleutels voor het losmaken van de deksel met schroefdraad die de holle doorgang van de motor bedekt (afbeelding 5.17). Schroef de deksel vervolgens los en verwijder deze.



Afbeelding 5.16. De veiligheidskap verwijderen.



Afbeelding 5.17. De dop met sleutels verwijderen.

- 3. Breng de aandrijfas in via de holle doorgang van de motoreenheid, zoals hierna beschreven. **Opmerking:** Smeer de uiteinden met schroefdraad van de uit meerdere segmenten bestaande aandrijfassen in met een lichte coating voor voeding goedgekeurde anti-seize bij elk gebruik. Wanneer u een uit meerdere segmenten bestaande aandrijfas installeert, moet u de segmenten met omgekeerde schroefdraad verbinden en de vergrendelpen van de deksel gebruiken om het segment tijdens het vastdraaien op zijn plek te houden.
 - Gebruik twee handen om de aandrijfas via de bovenkant van de mengeenheid te laden (afbeelding 5.18). Het inbrengen gaat gemakkelijker als u de as licht naar achteren en voren beweegt.
 - Wanneer er nog ongeveer 5,08 cm (2 inch) van de as resteert, draait u deze een beetje om de mengschroef vast te zetten.
 - Wanneer er nog ongeveer 2,54 cm (1 inch) van de as resteert, draait u deze een beetje om de lagereenheid vast te zetten.
 - Wanneer er nog ongeveer 0,64 cm (0,25 inch) van de as resteert, draait u deze om de spiebaan van de motoraandrijving uit te lijnen met de vier buitenste sleuven op de kop van de aandrijfas (afbeelding 5.19).





Afbeelding 5.18. Aandrijfas inbrengen. Afbeelding 5.19. Uitlijnen met de spiebaan.

- 4. Koppel de aandrijfas rechtstreeks aan de motoraandrijving.
 - Plaats de deksel met schroefdraad op de holle doorgang en draai deze met de hand rechtsom vast (afbeelding 5.20).
 - Plaats de moersleutel op de holle doorgang en draai de deksel vast met de bijgeleverde momentsleutel (afbeelding 5.21).
 Opmerking: De momentsleutel heeft een standaard vierkantaandrijving van 10 mm (3/8 inch) en is in de fabriek gekalibreerd op 16,9 Nm (150 in-lb.).
 - Controleer of de sleutels uit het systeem zijn verwijderd en in de opslaghouders zijn opgeborgen.
 - Sluit de veiligheidsdeksel. De magnetische sluiting houdt de veiligheidsdeksel gesloten.





Afbeelding 5.20. De dop terugplaatsen.

Afbeelding 5.21. De dop vastdraaien.

5.6 BPC's laden voor open-top mengen

Voor toepassingen met open-top mengen biedt de S.U.M. een herbruikbare lagerpoort (afbeelding 5.22), een mengschroef en mof voor eenmalig gebruik (afbeelding 5.23) en een tankvoering voor eenmalig gebruik met afvoerleiding.



Afbeelding 5.22. Lagerpotten voor drieklem (links) en snelaansluitingen (rechts) voor 2.000 l.





Voor open-top mengen in alle S.U.M.-tanks brengt u de herbruikbare lagerpoort (voor aansluiting met drie klemmen of snelaansluitingen) aan in de lagerpoortontvanger van de motor. De S.U.M. van 2.000 l gebruikt de versie met drie klemmen van de herbruikbare lagerpoort, terwijl alle andere maten de versie met snelaansluiting gebruiken (afbeelding 5.22). Er zijn ook twee passende mengschroef- en mofcombinaties: een met de aansluiting met drie klemmen en een met de snelaansluiting.

Volg de stappen onderstaande stappen om uw systeem voor opentop mengen te configureren.

 Laad de BPC-voering in de buitenste ondersteuningscontainer van de S.U.M. Volg de laadinstructies uit paragraaf 5.2 — De BPC laden, om de voering aan de hardware-eenheid te bevestigen. 2. Spreid de bovenste opening van de voering uit over de buitenrand van de hardware en hang de voering over de zijkanten (afbeelding 5.24).



Afbeelding 5.24. De voering installeren.

3. Open de houderklem voor de lager en breng de herbruikbare lagerpot in de ontvanger in (afbeelding 5.25). Sluit vervolgens de klem en vergrendel deze.



Afbeelding 5.25. De herbruikbare lagerpot inbrengen.

 Bevestig de mof van de mengschroef rechtstreeks op de herbruikbare lagerpoort met behulp van de snelaansluiting op de plastic huls. Een hoorbare klik geeft aan dat de bevestiging tot stand is gebracht (afbeelding 5.26). **Opmerking:** De eenheid van 2.000 I heeft geen snelaansluiting. Voor eenheden van 2.000 I wordt de verbinding tot stand gebracht met een aansluiting met drie klemmen.



Afbeelding 5.26. De mof van de mengschroef bevestigen.

5. Breng de aandrijfas aan in de motoras zoals beschreven in paragraaf 5.5 – De aandrijfas inbrengen.

5.7 Sondeaansluitingen maken

De instructies voor sondeaansluiting in deze paragraaf zijn van toepassing op Pall[™] Kleenpak[™] aseptische connectoren. **Opmerking:** Uw S.U.M.-systeem gebruikt in plaats van Kleenpak mogelijk CPC[™] AseptiQuik[™] aseptische connectoren of niet-aseptische snelaansluitingen. Neem contact op met uw salesmedewerker voor informatie over deze alternatieve sondeaansluitingen.

5.7.1 Kleenpak-specificaties

De Kleenpak-connector heeft een maximale werkdruk van 3 bar (43,5 psi) bij 40°C in compatibele vloeistoffen.



WAARSCHUWING: Een werking buiten de bovenstaande specificaties en/of met vloeistoffen die niet compatibel zijn met de bouwmaterialen kan leiden tot persoonlijk letsel of schade aan het apparaat.

5.7.2 Ontvangst van de uitrusting

De mannelijke en vrouwelijke Kleenpak-connectoren worden geleverd in afzonderlijke pakketten. Er zijn meerdere soorten eindaansluitingen om te voldoen aan de eisen van verschillende slangmaten en om verschillende aansluitmogelijkheden op slangen te bieden. Tips:

- Bewaar de mannelijke en vrouwelijke Kleenpak-connectoren in een schone, droge omgeving en zo mogelijk in de buitenverpakking waarin ze zijn afgeleverd.
- Verwijder de Kleenpak-connectoren NIET uit de binnenste apparaatzakverpakking waarin ze zijn geleverd.
- Mannelijke en vrouwelijke Kleenpak-connectoren worden beschermd geleverd in een binnen- en buitenzak. Controleer of de verpakking niet beschadigd is.
- De montagehulp wordt niet-steriel geleverd en kan meermaals worden gebruikt. De eenheid moet tussen elk gebruik worden bewaard in een schone, droge omgeving. Het montagehulpmiddel wordt apart geleverd en kan bij uw plaatselijke Pallvertegenwoordiger worden aangeschaft.

5.7.3 Installatie

Vóór installatie moet u absoluut controleren of de Kleenpak-connector geschikt is voor de vloeistof waarbij het voor de toepassing in aanraking komt, en moet u de betreffende onderstaande instructies volgen.

- Installeer de mannelijke en vrouwelijke Kleenpak-connectoren met compatibele verbindingen. Bevestig om lekken tijdens het gebruik te voorkomen de slang stevig aan de slangpilaar met kabelbinders of op een andere manier. Tijdens de montage van de buis wordt een vroegtijdige bediening van de mannelijke zuiger voorkomen door de tegenhoudring. De tegenhoudring moet op zijn plaats blijven totdat de feitelijke verbinding tot stand is gebracht. De aanwezigheid van kleppen vóór de connector op de slangen wordt aangeraden om te voorkomen dat vloeistof voorafgaand aan gebruik in aanraking komt met de Kleenpak-connectoren
- Als de connectoren moeten worden geautoclaveerd, plaatst u ze met de afpelstroken omhoog om blokkering ervan door condens te voorkomen.



WAARSCHUWING:

- Het apparaat moet droog blijven voor verbinding van de mannelijke en vrouwelijke Kleenpak-connectoren. Als er vloeistof in de leiding of rond de apparaten aanwezig is, mag u ze niet gebruiken.
- Deze Kleenpak-wegwerpconnectoren mogen niet met stoom gesteriliseerd worden in de leiding. Beperkingen inzake materiaalontwerp worden overschreden wanneer ze worden blootgesteld aan stoom onder druk en uiteindelijk zullen ze scheuren.

5.7.4 Gammastraling

- 1. Sluit de mannelijke of vrouwelijke Kleenpak-connector aan op het systeem voor eenmalig gebruik. Een klep of klem moet dicht bij de connector worden geïnstalleerd zodat ze niet nat worden nadat het systeem is gevuld met vloeistof.
- 2. Zorg dat de beschermkap goed is geplaatst. Autoclaafpapier en ander stralingsbestendig materiaal kan worden gebruikt om te vermijden dat de kap loskomt tijdens de hantering.
- 3. Het wordt aanbevolen om het volledige systeem in een binnen- en buitenzak te plaatsen ter bescherming vóór de gammastraling.
- 4. Behandel het met gammastraling. De maximaal toegelaten stralingsdosis is 50 kGy (5 mrad).

Belangrijke opmerking: Pall beveelt aan om de efficiëntie van de gammastralingscyclus te valideren met een geschikte methode. Deze connectoren zijn niet gevalideerd voor herhaalde blootstelling aan gammastraling.

5.7.5 Instructies voor het autoclaveren

- 1. Installeer de mannelijke of vrouwelijke Kleenpak-connector op de te autoclaveren uitrusting. Als de Kleenpak-connector op een tank is aangesloten, moet de tank juist worden geventileerd met een ontluchter.
- 2. Zorg dat de beschermkap van de Kleenpak-connector goed is geplaatst. Autoclaafpapier of een ander autoclaveerbaar en lucht/stoom doorlatend materiaal kan worden gebruikt om de dop losjes te bedekken zodat deze zich tijdens hantering niet kan verplaatsen.
- 3. De Kleenpak-connectoren moeten kunnen ventileren tijdens het autoclaveren. De ventilatiestrook moet naar boven wijzen om blokkering door condens te voorkomen.



WAARSCHUWING: Om ophoping van condens in de connectoren te voorkomen mag u de ventilatiestrook tijdens het autoclaveren niet naar beneden plaatsen. De connector mag niet worden afgedekt met zware voorwerpen tijdens de autoclaafcyclus. Pall beveelt aan om de efficiëntie van de autoclaveercyclus te valideren met een geschikte methode. **Opmerking:** De maximumtemperatuur is 121°C voor ACDonderdeelnummers en 130°C voor KPCHT-onderdeelnummers. De maximale blootstellingstijd is 75 minuten. Autoclaveer niet bij een hogere temperatuur of voor langere tijd. Een trage uitlaatcyclus wordt aanbevolen.

5.7.6 De verbinding tot stand brengen

Afbeelding 5.27 illustreert de onderdelen van een mannelijke en een vrouwelijke Kleenpak-connector en de volledige, geactiveerde connector.



WAARSCHUWING: Gebruik de connector niet als er vloeistof mee in aanraking komt. Gebruik de connector niet als de beschermdoppen los zitten of verplaatst zijn.



Afbeelding 5.27. Schema van Kleenpak-connector.

De verbinding tot stand brengen met het montagehulpmiddel voor connectoren

Volg de onderstaande stappen om de Kleenpak-verbinding tot stand te brengen met het montagehulpmiddel. 1. Trek het lipje omhoog en verwijder deze van de beschermdoppen zodat u de doppen kunt verwijderen van de Kleenpakconnectoren (afbeelding 5.28).



Afbeelding 5.28. Beschermdoppen op connectoren.

 Houd de cilinder van de grote (mannelijke) Kleenpak-connector boven de basis. Lijn de kleine (vrouwelijke) Kleenpak-connector uit met de mannelijke connector. De platte zijden moeten worden uitgelijnd en beide afpelbare stroken moeten gevouwen blijven (afbeelding 5.29). **Opmerking:** Indien de Kleenpakconnectoren niet juist zijn uitgelijnd, kan geen verbinding tot stand worden gebracht.



Afbeelding 5.29. Uitgelijnde platte kanten van connectoren.

3. Nadat beide connectoren juist zijn uitgelijnd, drukt u ze stevig tegen elkaar totdat beide vergrendelclips stevig dichtklikken (afbeelding 5.30).



Afbeelding 5.30. Connectoren samendrukken.

 Ondersteun de mannelijke en vrouwelijke Kleenpak-connectoren en verwijder de tegenhoudring uit de mannelijke connector door het lipje naar het getande uiteinde van de mannelijke Kleenpakconnector te trekken (afbeelding 5.31).



Afbeelding 5.31. De tegenhoudring verwijderen.

5. Als u het montagehulpmiddel voor Kleenpak-connectoren gebruikt, plaatst u de connector zodanig in het hulpmiddel dat de afpelbare stroken door de opening naar buiten steken (afbeelding 5.32). Als dat niet zo is, houdt u de connector stevig vast aan beide zijden, waarbij de afpelbare stroken van uw hand af wijzen. De Kleenpak-connector moet goed in het montagehulpmiddel blijven zitten als deze juist is geïnstalleerd.





6. Houd het montagehulpmiddel in de palm van uw hand. Hierbij moet de Kleenpak-connector naar buiten wijzen en moet u met uw duim de Kleenpak-connector in het montagehulpmiddel ondersteunen. Pak met uw andere hand beide afpelbare stroken zo dicht mogelijk bij de behuizing van het montagehulpmiddel vast om een goede grip te garanderen, en trek beide stroken in één vloeiende beweging los. Zorg dat de Kleenpak-connector loodrecht op de afpelbare stroken staat (afbeelding 5.33).



Afbeelding 5.33. De afpelbare stroken van de connector verwijderen.



WAARSCHUWING: Gebruik de connector niet als maar één afpelbare strook is verwijderd in plaats van beide.

 Terwijl de Kleenpak-connector nog stevig in het montagehulpmiddel zit, duwt u de duimsteun van de mannelijke Kleenpak-connector naar de basis van de cilinder (afbeelding 5.34).



Afbeelding 5.34. De duimsteun van de connector naar de basis van de cilinder duwen.

Voor een goede verbinding moet de plunjer in de mannelijke Kleenpak-connector volledig in de vrouwelijke Kleenpak-connector zijn ingebracht. Bij wijze van controle herhaalt u de activering totdat u een harde stop bereikt.

U kunt de Kleenpak-connector zo nodig uit het montagehulpmiddel verwijderen om de plunjerbeweging te voltooien. 8. Na assemblage van de Kleenpak-connector kan de montagehulp worden verwijderd. Wanneer de montagehulp wordt verwijderd, verifieert u de bediening tot een harde aanslag is bereikt. Start vervolgens de vloeistofoverdracht (afbeelding 5.35)



Afbeelding 5.35. Vloeistofoverdracht beginnen.

De verbinding tot stand brengen zonder het montagehulpmiddel voor connectoren

Volg de onderstaande stappen om de Kleenpak-verbinding tot stand te brengen zonder het montagehulpmiddel.

1. Trek de beschermdoppen omhoog zodat u de doppen kunt verwijderen van de Kleenpak-connectoren (afbeelding 5.36).



Afbeelding 5.36. Beschermdoppen verwijderen.

 Houd de cilinder van de grootste (mannelijke) connector boven de basis. Lijn de kleinste (vrouwelijke) connector uit met de mannelijke connector. De platte zijden moeten worden uitgelijnd en beide afpelbare stroken moeten gevouwen blijven (afbeelding 5.37).
 Opmerking: Indien de Kleenpak-connectoren niet juist zijn uitgelijnd, kan geen verbinding tot stand worden gebracht.



Afbeelding 5.37. Mannelijke en vrouwelijke connectoren uitlijnen.

3. Nadat beide connectoren juist zijn uitgelijnd, drukt u ze stevig tegen elkaar totdat beide vergrendelclips stevig dichtklikken (afbeelding 5.38).



Afbeelding 5.38. De connectoren samendrukken.

4. Ondersteun zowel de mannelijke en als de vrouwelijke Kleenpakconnector. Verwijder de tegenhoudring van de mannelijke connector door de tab naar het geribbelde uiteinde van de mannelijke Kleenpak-connector te trekken (afbeelding 5.39).



Afbeelding 5.39. De tegenhoudring verwijderen.

5. Ondersteun met één hand de mannelijke en vrouwelijke kanten van de Kleenpak-connector door uw vingers rond beide kanten van de connector te wikkelen, naast de flens. Pak met uw andere hand beide afpelbare stroken zo dicht mogelijk bij de platte kant van de connector vast om een goede grip te garanderen, en trek ze in één vloeiende beweging uit elkaar. Zorg dat de connector loodrecht op de trekstrips staat, zoals getoond in afbeelding 5.40. De loodrechte oriëntatie moet aangehouden worden wanneer tegelijk aan de twee strips wordt getrokken. Niet gebruiken indien een van de trekstrips is verwijderd in plaats van beide.

Opmerking: Oefen geen loodrechte druk uit op de connector omdat die dan kan breken. Als er een loodrechte kracht op wordt uitgeoefend door items die aan de Kleenpak-connector zijn bevestigd, moet de connector juist worden ondersteund.



Afbeelding 5.40. De afpelbare stroken van de connector verwijderen.

6. Duw de duimsteun van de mannelijke Kleenpak-connector naar de basis van de cilinder totdat ze elkaar raken (afbeelding 5.41).



Afbeelding 5.41. De duimsteun naar de cilinder duwen.

Voor een goede verbinding moet de plunjer in de mannelijke connector volledig in de vrouwelijke connector zijn ingebracht. Bij wijze van controle herhaalt u de activering totdat u een harde stop bereikt. Start vervolgens de vloeistofoverdracht (afbeelding 5.42)



Afbeelding 5.42. Vloeistofoverdracht beginnen.

5.8 Sensoren en sondes inbrengen

5.8.1 Temperatuursensoren inbrengen

Volg de onderstaande stappen om temperatuursensoren in de thermowell in te brengen.

- 1. Verwijder het plastic inzetstuk in de dompelhuls, indien aanwezig.
- Plaats een kleine hoeveelheid glycerol (0,5 ml) in de dompelhuls om te helpen bij de warmteoverdracht. De glycerol dient ook als smeermiddel en helpt bij het inbrengen. Draai de temperatuursensor lichtjes terwijl u hem begint in te brengen (afbeelding 5.43).



Afbeelding 5.43. Temperatuursensor inbrengen.

3. Breng de temperatuursensor in totdat de basis van de sonde de mond van de thermowell raakt.

4. Zet de sensor vast door de kraag van de Luer-lock te verdraaien (afbeelding 5.44). De thermowell rekt iets op wanneer de sensor wordt aangebracht.



Afbeelding 5.44. De temperatuursensor verdraaien om deze goed aan te brengen.

5.8.2 pH- en geleidbaarheidssondes voorbereiden

- Selecteer de gewenste sonde. Zorg dat er een Teflon[™]ondersteuningsring en een O-ring op de sonde aanwezig zijn en
 inspecteer ze op het oog op beschadiging.
- Voer eventueel vereist onderhoud van de sondes uit en kalibreer de pH- of geleidbaarheidssonde (hoofdstuk 4 – Kalibratieprocedures).
- 3. Doe een klein beetje (0,5–1 ml) buffer- of fysiologische zoutoplossing op het uiteinde met schroefdraad van de sondeadapter.
- 4. Steek de sonde in de sonde-eenheid door de adapter met schroefdraad.
- Zorg dat de sondetip het membraan van de aseptische connector of de snelaansluiting niet raakt door een ruimte van meer dan 6,35 mm (1/4 inch) vrij te laten. Schroef de sondetip vervolgens in de sondeadapter.
- 6. Draai de adapter met de hand vast en verifieer dat de sondepunt het membraan niet raakt.
- Plaats de sonde-eenheid met de sonde in de autoclaaftray voor sondekits (afbeelding 5.45). Opmerking: Afbeelding 5.45 laat zien hoe de autoclaaftray wordt gebruikt voor sondes met Kleenpak connectoren. Voor sondes met AseptiQuik aseptische connectoren is een andere autoclaaftray vereist.



Afbeelding 5.45. Kleenpak-sondeeenheid in autoclaaftray.

- Autoclaveer de sonde-eenheid volgens een gevalideerde sterilisatiecyclus (ongeveer 30 minuten op 122°C). Een sterilisatiecyclus van 30 minuten is doorgaans voldoende. Opties voor droge of natte parameters kunnen worden gebruikt. Trage uitlaatcycli hebben de voorkeur, aangezien dit de sondes minder belast tijdens temperatuur- en drukwijzigingen van het autoclaveren.
- Indien juist bewaard kunnen de geautoclaveerde sondeeenheden kortstondig (minder dan 24 uur) droog worden bewaard zonder dat de levensduur, prestaties of steriliteit van de sensor achteruit gaat.

Opmerking: Voordat u een sonde inbrengt, dient u zich vertrouwd te maken met de instructies voor Kleenpak-connectoren in paragraaf 5.7.

5.8.3 pH- en geleidbaarheidssondes inbrengen

Volg de onderstaande stappen om de pH- en geleidbaarheidssondes in de BPC in te brengen met behulp van de sonde-eenheid (geautoclaveerd indien vereist). **Opmerking:** Afbeeldingen 5.47 en 5.48 tonen een plastic sondeclip. Uw systeem is mogelijk al uitgerust met een metalen sondeclip.

 Plaats slangklemmen op alle sondepoorten voordat u probeert sonde-eenheden aan te sluiten (afbeelding 5.46). Dit voorkomt verlies van steriliteit als er tijdens de aseptische aansluiting fouten worden gemaakt.



Afbeelding 5.46. Slangklemmen plaatsen.

- 2. Sluit de aseptische connector aan op de poort voor pH- of geleidbaarheidssondes. **Opmerking:** Als u Kleenpak aseptische connectoren gebruikt, volgt u de methodes uitgelegd in paragraaf 5.7.
- 3. Zorg dat alle vier de klemmen vastklikken op elke connector en dat de basis en cilinder volledig zijn ingebracht.
- 4. Verwijder de slangklem op de individuele sondepoort.
- Breng de pH/geleidbaarheidssonde in door de balgen naar de S.U.M. samen te vouwen. (afbeelding 5.47). De sonde is goed ingebracht wanneer de balgen volledig zijn samengevouwen. Een klein beetje vloeistof kan in de balg binnendringen als de sonde wordt geladen na het vullen met vloeistof.



Afbeelding 5.47. De sondebalgen samendrukken.

- 6. Bevestig een sondeclip op de horizontale band onderaan op de voorkant van de buitenste ondersteuningscontainer om de sonde te ondersteunen.
- 7. Plaats de volledig samengevouwen sonde-eenheid in de sondeclip (afbeelding 5.48). Druk de sondebalgen zo nodig verder samen om de sonde-eenheid stevig in het uiteinde van de sondeclip aan te brengen.



Afbeelding 5.48. De samengevouwen sonde-eenheid in de sondeclip plaatsen.

8. Laat de sondebalgen ontspannen.

Opmerking: Sondepoorten die niet worden gebruikt voordat vloeistof wordt toegediend, moeten afgeklemd blijven. Niet-gebruikte sondepoorten zijn later toegankelijk om sondes te vervangen of voor gebruik als redundante poort, mocht dat nodig zijn.

5.9 De S.U.M. tarreren

Om nauwkeurige massametingen te waarborgen moet u uw S.U.M.systeem met behulp van de volgende stappen tarreren voordat u de BPC vult met vloeistof. **Opmerking:** Als u laadcellen gebruikt, moet u de module Mass (Massa) op de touchscreenconsole configureren voordat u het systeem kunt tarreren. Zie paragraaf 3.5.3 voor informatie over het configureren van de module Mass (Massa).

 Controleer of alle onderdelen van de mixer (aandrijfas, BPC, sondedeurplaten en gereedschappen) zich op de buitenste ondersteuningscontainer bevinden. De BPC moet ook in de eenheid zijn aangebracht. Afhankelijk van uw weegprocessen moet de Powdertainer mogelijk ook worden geïnstalleerd.

- 2. als u geen laadcellen gebruikt, tarreert u de skid op de schaal terwijl de BPC is geïnstalleerd.
- Als u laadcellen gebruikt, moet de module Mass (Massa) op de touchscreenconsole zijn geconfigureerd. Raak op het startscherm de module Mass (Massa) aan en selecteer 'Calibration' (Kalibratie). Raak '1-Point Calibration' (1-punts kalibratie) aan. Voer 0 (nul) in het veld in en raak 'Next' (Volgende) aan. De S.U.M. is getarreerd wanneer het scherm '1 Point Calibration Complete' (1-punts kalibratie voltooid) verschijnt.

6

Bedieningsinformatie

Inhoudsopgave

6.1

- Vullen met vloeistof
- 6.2 Roeren instellen
- 6.3 Poeder toevoegen met de Powdertainer
- 6.4 De timer gebruiken
- 6.5 Functies voor bewaking en controle
- 6.6 Aftappen
- 6.7 Voorbereiden op de volgende run

De instructies in dit hoofdstuk zijn een algemeen overzicht van de bedieningsprocedures. Omdat procedures per toepassing sterk kunnen verschillen, is het belangrijk uw interne formuleringsprotocollen te raadplegen voor de juiste procedures.

De onderstaande informatie heeft betrekking op steriele en nietsteriele mengtoepassingen. Voor steriele toepassingen moeten alle aansluitingen aseptisch worden gemaakt.

6.1 Vullen met vloeistof

Als uw systeem laadcellen en pompen heeft die zijn geconfigureerd zoals in paragraaf 3.1, volgt u de onderstaande stappen om het systeem te vullen met de module Fill (Vullen) op de touchscreenconsole. **Opmerking:** De module Fill (Vullen) moet vóór gebruik zijn geconfigureerd. Zie paragraaf 3.5.7 voor informatie over het configureren van de module Fill (Vullen).

- Zoek de module Fill (Vullen) op het startscherm van de touchscreenconsole op. Raak de module aan om het scherm Fill Settings (Vulinstellingen) te openen en selecteer vervolgens de knop 'Fill control' (Vullen reguleren).
- 2. In het scherm Fill control (Vullen reguleren) wordt u gevraagd een vulmethode te selecteren: door middel van druk of een pomp.
 - Als u pressure (druk) selecteert: Voer het vloeistofdoel in. Raak vervolgens 'Done' (Gereed) aan.
 - Als u pump (pomp) selecteert: Voer het vloeistofdoel in door de hoeveelheid in te voeren die aan de huidige waarde moet worden toegevoegd. Voer de stroomsnelheid voor fase 1 in, het percentage gevuld vóór fase 2 en de stroomsnelheid voor fase 2. Raak vervolgens 'Done' (Gereed) aan.
- 3. Als het systeem aangesloten knijpventielen bevat, wordt u in het volgende geopende scherm gevraagd te controleren of het knijpventiel voor vullen is gesloten. Als het knijpventiel open is, verplaatst u de schuifknop van 'Open' naar 'Closed' (Gesloten). Raak vervolgens 'Start Fill' (Vullen starten) aan.
- 4. Nadat u 'Start Fill' (Vullen starten) hebt geselecteerd, worden de pompen ingeschakeld en begint het vullen van het systeem met de door de gebruiker gespecificeerde hoeveelheid vloeistof. Het geopende scherm (afbeelding 6.1) toont de resterende hoeveelheid vloeistof, de huidige massa, de BPC-druk en de vloeistofdruk (voor vloeistofdruk1, vloeistofdruk2 en delta).

 U kunt te allen tijde onder aan het scherm de knop 'Pause' (Pauzeren) selecteren om het vulproces tijdelijk te stoppen. Er verschijnt een scherm Fill Paused (Vullen gepauzeerd) waarmee u het vullen met vloeistof kunt hervatten of annuleren.



Afbeelding 6.1. Het scherm Fill Control (Vullen reguleren).

6.2 Roeren instellen

Wanneer de vloeistof het minimale gebruiksvolume van de mixer voor eenmalig gebruik (S.U.M.) heeft bereikt, kunt u het roeren inschakelen met de module Agitation (Roeren) op de touchscreenconsole. Zorg dat de BioProcess Container (BPC) tot een adequaat volumeniveau voor uw individuele lopende processen is gevuld voordat u met roeren begint.

VOORZICHTIG: Als u de S.U.M. buiten de in dit document vermelde mengspecificaties gebruikt, kan de aandrijfas beschadigd raken.

Opmerking: De module Agitation (Roeren) moet zijn geconfigureerd voordat u begint met mengen. Zie paragraaf 3.5.2 voor informatie over het configureren van de module Agitation (Roeren). Als u het roeren wilt inschakelen, zoekt u de module Agitation (Roeren) op het startscherm van de touchscreenconsole op (afbeelding 6.2) en raakt u de knop 'Start' (Starten) aan.



Afbeelding 6.2. Module Agitation (Roeren).

6.3 Poeder toevoegen met de Powdertainer

De S.U.M. kan worden uitgerust met een Powdertainerondersteuningsarm om vaste stoffen zoals poeder toe te voegen. Zie paragraaf 8.5.2 van deze publicatie voor specificaties van de Powdertainer.

In de volgende instructies wordt het gebruik van de Powdertainer-BPC in combinatie met de S.U.M. beschreven wanneer deze is voorzien van de Powdertainer-arm. Hoewel er ook andere methodes zijn om vaste stoffen met de BPC-poederpoort toe te voegen, zijn hier geen volledige protocollen opgenomen omdat individuele protocollen kunnen wisselen.

- 1. Zorg dat de klem op de zak van de Powdertainer volledig gesloten is.
- 2. Hang de Powdertainer met de handgreep aan de hanger van de ondersteuningsarm (afbeelding 6.3).



Afbeelding 6.3. De Powdertainer aan de ondersteuningsarm hangen.

- 3. Pas de hoogte van de Powdertainer-arm zo nodig aan door de vergrendeling omlaag te klappen, de bovenste stang omhoog of omlaag te schuiven, de vergrendeling opnieuw te sluiten en de hendel vervolgens rechtsom te draaien om deze vast te zetten.
- 4. Verwijder de tape en dop van de opening van de Powdertainer-BPC (afbeelding 6.4).



Afbeelding 6.4. De dop van de Powdertainer verwijderen.

- Klik de poederpoort op de BPC in de beugel op de ondersteuningsarm van de Powdertainer. Verplaats de beugel zo nodig naar binnen of buiten om deze aan de positie van de poederpoort op de BPC aan te passen.
- 6. Verwijder de klem en beschermdop van de poederpoort op de BPC.
- 7. Lijn de opening van de Powdertainer en de poederpoort op de BPC met elkaar uit (afbeelding 6.5).



Afbeelding 6.5. BPC-poederpoort in de klem.

8. Gebruik de aansluiting met drie klemmen op de Powdertainer-arm om de twee poorten te verbinden (afbeelding 6.6).



Afbeelding 6.6. De Powdertainer aansluiten op de BPC-poederpoort.

9. Als u wilt beginnen met het toevoegen van vaste stoffen, maakt u de klem op de onderkant van de toevoerzak los (afbeelding 6.7). Zorg dat de volledige inhoud van de toevoerzak is toegevoegd voordat u de Powdertainer losmaakt van de BPC. De afdekking, pakking en klem van de poederpoort kunnen worden teruggeplaatst als er geen andere droge ingrediënten hoeven te worden toegevoegd, of de poederpoort kan open blijven voor monstername met een pipet.



Afbeelding 6.7. De Powdertainer-toevoerklem openen.

6.4 De timer gebruiken

Met de module Timer op de touchscreenconsole kunnen gebruikers een timer instellen om te bewaken hoe lang de S.U.M. aan het mengen is geweest. Zoek de module Timer op het startscherm van de touchscreenconsole op om de timer te gebruiken. Raak de knop 'Start' (Starten) aan. De timer begint te tellen vanaf nul en gaat door totdat de gebruiker de knop 'Stop' (Stoppen) aanraakt.

6.5 Functies voor bewaking en controle

6.5.1 Mengen

De gebruiker kan de mengsnelheid tijdens het gebruik bewaken met de module Agitation (Roeren) op de touchscreenconsole. De module Agitation (Roeren) op het startscherm toont de huidige mengsnelheid en de instelpunten die tijdens het configureren van de module zijn opgegeven. Als u de roersnelheid wilt kalibreren, raadpleegt u paragraaf 4.4 van deze publicatie.
6.5.2 Meer poeder of vloeistof toevoegen

Afhankelijk van uw bedieningsprocedures moet u tijdens het mengproces mogelijk meer poeder of vloeistof aan de oplossing toevoegen. Volg de onderstaande stappen om tijdens het gebruik componenten toe te voegen aan de S.U.M.

Poeder toevoegen aan vloeistof

- 1. Hang de Powdertainer-BPC aan de hanger op de S.U.M.
- 2. Verwijder de klem, afdekking en pakking van de poederpoort op de BPC.
- 3. Gebruik de houderarm om de poederpoort van de BPC op zijn plek te houden terwijl de Powdertainer wordt gepositioneerd.
- 4. Verwijder de beschermtape en -dop van de poort op de Powdertainer.
- 5. Sluit de Powdertainer met de aansluiting met drie klemmen aan op de poederpoort op de BPC.
- Controleer in de module Agitation (Roeren) op de touchscreenconsole of de motor van de mixer werkt op de gewenste snelheid. Zie paragraaf 6.5.1 voor meer informatie over het gebruik van de module Agitation (Roeren).
- 7. Open de clip van de zak om de inhoud van de Powdertainer in de S.U.M. BPC te laten lopen.
- 8. Blijf mengen totdat alle poeder is opgelost en voordat u doorgaat naar het volgende ingrediënt.

Vloeistof zonder pompen toevoegen aan vloeistof

- Controleer in de module Agitation (Roeren) op de touchscreenconsole of de motor van de mixer werkt op de gewenste snelheid. Zie paragraaf 6.5.1 voor meer informatie over het gebruik van de module Agitation (Roeren).
- 2. Sluit de leiding van de vloeistofbron met de snelaansluiting of aansluiting met drie klemmen aan op een extra leidingenset op de bovenkant van de BPC.
- Om de vloeistof toe te dienen opent u het vloeistoftraject door de klemmen op de leidingenset te openen. Blijf mengen totdat alle vloeistof is toegediend en voordat u doorgaat naar het volgende ingrediënt.

Vloeistof met pompen toevoegen aan vloeistof

- Controleer in de module Agitation (Roeren) op de touchscreenconsole of de motor van de mixer werkt op de gewenste snelheid. Zie paragraaf 6.5.1 voor meer informatie over het gebruik van de module Agitation (Roeren).
- Zorg dat de gebruikte pomp is geconfigureerd volgens de instructies in paragraaf 3.1 en dat de juiste pomp is toegewezen aan de module Fill (Vullen). Raak de module Fill (Vullen) aan op het startscherm van de touchscreenconsole om het scherm Fill Settings (Vulinstellingen) te openen. Raak de knop 'Pump Selection' (Pomp selecteren) aan om de pomptoewijzingen te controleren en/of bewerken.
- Ga terug naar het scherm Fill Settings (Vulinstellingen). Raak de knop 'Fill control' (Vullen reguleren) aan om de hoeveelheid vloeistof die u wilt toedienen, de fases, de stroomsnelheden en de overgangs- en beëindigingsparameters op te geven. Raak de knop 'Done' (Gereed) aan.
- 4. In het volgende scherm wordt u gevraagd of u knijpventielen gebruikt.
 - Als dat niet zo is: Raak de knop 'No' (Nee) aan. U wordt in een scherm gevraagd of u wilt beginnen met het vullen met vloeistof. Raak de knop 'Cancel' (Annuleren) aan om terug te gaan naar het vorige scherm, of raak 'Start Fill' (Vullen starten) aan om met vullen te beginnen.
 - Als dat zo is: Raak de knop 'Yes' (Ja) aan. Er verschijnt een scherm om de knijpventielen te openen of sluiten. Het knijpventiel moet worden gesloten om te beginnen met het vullen met vloeistof. Raak de knop 'Cancel' (Annuleren) aan om terug te gaan naar het vorige scherm, of raak 'Start Fill' (Vullen starten) aan om met vullen te beginnen.
- Nadat u 'Start Fill' (Vullen starten) hebt geselecteerd, worden de pompen ingeschakeld en begint het vullen van het systeem met de door de gebruiker gespecificeerde hoeveelheid vloeistof. Het geopende scherm toont de vloeistofdruk en deltawaarden, de resterende hoeveelheid vloeistof, de huidige massa en de BPC-druk.
- U kunt te allen tijde onder aan het scherm de knop 'Pause Fill' (Vullen pauzeren) selecteren om het vulproces tijdelijk te stoppen. Er verschijnt een scherm Fill Paused (Vullen gepauzeerd) waarmee u het vullen met vloeistof kunt hervatten of annuleren.

7. Ga door met mengen totdat alle vloeistof is toegediend. Ga daarna verder met het volgende ingrediënt, indien nodig.

Herhaal dit proces voor alle vloeistofcomponenten in de formulering.

6.5.3 pH en geleidbaarheid aanpassen

Op de touchscreenconsole kunnen gebruikers de pH- en geleidbaarheidsniveaus van de oplossing tijdens het mengen bewaken. Op het startscherm van de touchscreenconsole tonen de modules pH en Conductivity (Geleidbaarheid) de huidige waarden die door het systeem worden afgelezen. Als u deze niveaus wilt bijstellen, moet u de pompen gebruiken die zijn geconfigureerd voor zuur-, base- of fysiologische zoutoplossingen. Volg de onderstaande stappen om de pH- en geleidbaarheidsniveaus aan te passen.

- Voordat u de pH of geleidbaarheid aanpast, moet u zorgen dat de door u gebruikte sonde juist is gekalibreerd en goed is ingebracht in de sondebalgen. Zie paragrafen 4.2 en 4.3 voor informatie over de kalibratie van pH en geleidbaarheid.
- 2. Raak op het startscherm van de touchscreenconsole de module 'Pump' (Pomp) aan die is toegewezen aan de buffer die u wilt toedienen, zoals zuur, base of fysiologische zoutoplossing.
- 3. Selecteer 'Add Bolus' (Bolus toevoegen). In het geopende scherm wordt u gevraagd de gewenste eenheden, de hoeveelheid vloeistof die de pomp moet toedienen, en de toedieningsduur (duur van bolustoediening) op te geven. Het veld Set point flow rate (Stroomsnelheid instelpunt) wordt automatisch ingevuld nadat u de vloeistoftoediening en de duur hebt opgegeven. Opmerking: Als een duur of vloeistof wordt opgegeven die groter is dan de waarden waarvoor de pomp is gekalibreerd, wordt in deze velden automatisch de maximaal mogelijke waarde ingevuld. Als u geen controlevergelijking gebruikt, raakt u 'Next' (Volgende) aan om door te gaan. Als u parameters voor pH of geleidbaarheid wilt opgeven met een controlevergelijking, volgt u de onderstaande stappen voordat u de knop 'Next' (Volgende) aanraakt.
 - Schuif de aan/uitknop voor 'Control Equation' (Controlevergelijking) en 'Hysteresis' (Hysterese) naar 'On' (Aan).
 - Raak het veld naast 'Control Equation' (Controlevergelijking) aan om de instelpunten voor pH of geleidbaarheid op te geven en om op te geven welke actie moet worden ondernomen (Pause (Pauzeren) of Stop (Stoppen)).

- Raak het veld naast 'Hysteresis' (Hysterese) aan om de te bereiken pH- of geleidbaarheidswaarde op te geven en hoe lang die waarde moet worden aangehouden voordat de pomp doorgaat. Raak 'Next' (Volgende) aan.
- 4. Bovenaan in het geopende scherm wordt het volgende weergegeven: de resterende hoeveelheid toe te dienen vloeistof, de huidige pH- of geleidbaarheidswaarde en de totale hoeveelheid toe te dienen vloeistof. De resterende tijd is zichtbaar in het midden van het scherm. Als u een controlevergelijking hebt opgegeven, wordt de vergelijking onder aan het scherm weergegeven. Als er geen controlevergelijking is opgegeven, wordt in plaats daarvan onder aan het scherm 'Control Equation Disabled' (Controlevergelijking uitgeschakeld) weergegeven.

6.5.4 Temperatuur bijstellen

Op de touchscreenconsole kunnen gebruikers de temperatuur van de oplossing tijdens het mengen bewaken. Op het startscherm van de touchscreenconsole toont de module Temperature (Temperatuur) de huidige temperatuur in graden Celsius (°C) die door het systeem wordt afgelezen. Als u de temperatuur van de S.U.M. wilt instellen met de module Temperature (Temperatuur) op de touchscreenconsole, moet op uw systeem een geïntegreerde temperatuurcontrole-unit (TCU) zijn geconfigureerd. Volg de onderstaande instructies om de S.U.M.-temperatuur aan te passen.

- 1. Voordat u de temperatuur aanpast, moet u zorgen dat de door u gebruikte sonde juist is ingebracht in de sondebalgen.
- Raak op het startscherm van de touchscreenconsole de module Temperature (Temperatuur) aan. Raak de knop 'Select Active Sensor' (Actieve sensor selecteren) aan op het scherm Temperature Settings (Temperatuurinstellingen).
- In het volgende scherm wordt de gebruiker gevraagd een temperatuursensor te selecteren. U kunt Temp1 of Temp2 kiezen. Nadat u een sensor hebt geselecteerd, raakt u 'Done' (Gereed) aan. Raak 'Close' (Sluiten) aan om terug te gaan naar het startscherm van de touchscreenconsole.

4. Raak de knop 'Setpoint' (Instelpunt) op de module Temperature (Temperatuur) aan. Geef het gewenste TCU-instelpunt (in°C) op en raak 'Enter' aan. Raak 'Start' (Starten) op de module Temperature (Temperatuur) aan. De TCU past nu de temperatuur aan op grond van het verschafte instelpunt. De PID-luswaarden voor verwarmen en koelen zijn in de TCU opgenomen.

6.5.5 De druk in de BPC-hoofdruimte bewaken

Op de touchscreenconsole kunnen gebruikers de druk in de hoofdruimte van de BPC bewaken tijdens het gebruik van de S.U.M. De module BPC Pressure (BPC-druk) op het startscherm van de touchscreenconsole (afbeelding 6.8) toont de huidige BPC-druk in psi.



Afbeelding 6.8. BPCdrukmodule.

6.5.6 Alarmen bewaken

In veel van de modules op de touchscreenconsole kunnen gebruikers de alarmen High/Low (Hoog/Laag) of High High (Hoog-hoog)/Low Low (Laag-laag) instellen. Deze alarmen worden geactiveerd onder door de gebruiker gespecificeerde voorwaarden. Deze voorwaarden worden binnen individuele modules ingesteld.

Alarmen High (Hoog) en Low (Laag)

Als een alarm High (Hoog) of Low (Laag) tijdens het gebruik wordt geactiveerd:

- De betreffende module wordt geel op het startscherm.
- De gebruikers worden visueel gewaarschuwd voor specifieke omstandigheden, maar het functioneren wordt niet gepauzeerd of gestopt.

Alarmen High High (Hoog-hoog) en Low Low (Laag-laag)

Als een alarm High High (Hoog-hoog) of Low Low (Laag-laag) tijdens het gebruik wordt geactiveerd:

- De betreffende module wordt rood op het startscherm.
- Er klinkt een hoorbare zoemer.
- Operationele processen die zijn gerelateerd aan de betreffende module worden gepauzeerd of gestopt, afhankelijk van de door de gebruiker gespecificeerde voorwaarden.

Geactiveerde alarmen deactiveren

Als u een geactiveerd alarm wilt deactiveren, raakt u het alarmpictogram aan op het startscherm van de touchscreenconsole om het scherm Alarm Status (Alarmstatus) te openen. Raak het geactiveerde alarm aan (of selecteer de knop 'Select All' (Alles selecteren) als u meerdere alarmen tegelijk wilt bevestigen) en raak vervolgens 'Acknowledge' (Bevestigen) aan. De zoemer stopt en het alarmpictogram van de betreffende module verandert van rood terug in wit.

6.5.7 Andere functies bewaken

U kunt de touchscreenconsole gebruiken om andere functies buiten de geleverde modules te bewaken. Met de hierna beschreven modules Auxiliary Output (Hulpuitgang) en Auxiliary Input (Hulpingang) kunnen gebruikers aanpassen welke functies ze willen bewaken tijdens het mengproces. Beheerders hebben bovendien toegang tot een optie 'Diagnostics' (Diagnose) in de instellingen voor de pomp- en pH-modules.

Hulpuitgangen

In de module Auxiliary Output (Hulpuitgang) kunnen gebruikers een bestaande module dupliceren om over extra sensoren of meer redundantie te beschikken. Als u geconfigureerde hulpuitgangen tijdens het gebruik wilt bewaken, raakt u op het startscherm van de touchscreenconsole de module Auxiliary Output (Hulpuitgang) aan. Nadat u een module hebt geselecteerd die u wilt dupliceren, kunt u een diagnose kiezen, de modulenaam wijzigen en de RPM-schaling aanpassen.

Hulpingangen

In de module Auxiliary Input (Hulpingang) kunnen gebruikers sensoren en zenders bewaken die een signaal van 4-20 mA gebruiken en geen gespecificeerde module op de touchscreenconsole hebben.

Pomp- en pH-diagnose

Beheerders hebben toegang tot de optie 'Diagnostics' (Diagnose) voor de pomp- en pH-modules. Met Pump Diagnostics (Pompdiagnose) kan de beheerder de uitgangsstroom van de Pompuitgang handmatig instellen. Dit kan helpen de stroom naar RPM-waarden voor RPM-schaling te bepalen. Met pH Diagnostics (pH-diagnose) kan de beheerder de kalibratiewaarden voor de helling en offset weergeven die momenteel op de pH-sensor(en) worden toegepast.

6.5.8 Monstername

Monstername met recirculatielus

Monsters kunnen worden genomen uit de SmartSite[™]-poort op de recirculatielus. U brengt daarvoor een injectiespuit met Luer-lock aan en neemt een monster af terwijl de recirculatielus in gebruik is. U kunt ook monsters nemen op BPC's die zijn uitgerust met een thermowell-monsternamepoort in een van de sondepoorten. De volgende instructies zijn bedoeld voor aseptische monstername via de SmartSite-poort. Gebruik voor deze instructies de standaard Luerlock op een injectiespuit van 60 ml of een steriel verdeelstuk.

- 1. Verwijder het stof van de SmartSite (afbeelding 6.9).
- 2. Reinig de SmartSite met een nat doekje. Sluit vervolgens de sanitaire injectiespuit met Luer-Lock aan (afbeelding 6.10).





Afbeelding 6.9. Dop verwijderen.

Afbeelding 6.10. Injectiespuit voor monstername inbrengen.

- Pas een beetje vacuümdruk toe door de plunjer van de injectiespuit iets naar buiten te trekken en een monster te nemen (circa 30-60 ml)
- 4. Verwijder de injectiespuit. Dit wordt gezien als een zuiveringsmonster.
- 5. Reinig de SmartSite met een nat doekje. Sluit vervolgens de sanitaire injectiespuit met Luer-Lock aan.
- 6. Schep een lichte vacuümdruk met de spuit en neem het gewenste monster (ongeveer 10–20 ml).
- 7. Verwijder de injectiespuit. Dit wordt gezien als een representatief monster.
- 8. Reinig de SmartSite met een nat doekje. Plaats de stofdop vervolgens weer terug.

Monstername zonder recirculatielus

Wanneer de recirculatielus niet wordt gebruikt, kan in plaats daarvan een leidingenset worden gebruikt voor monstername. Als u een leidingenset voor monstername wilt gebruiken, duwt u de leidingenset omlaag tot het vloeistofniveau. Gebruik een pomp om vloeistof af te tappen naar een secundaire container. Monsters kunnen op dezelfde manier ook worden genomen via de afvoerleiding.

6.6 Aftappen

6.6.1 De module Harvest (Aftappen) gebruiken

Als u pompen hebt geconfigureerd zoals in paragraaf 3.1, volgt u de onderstaande stappen om het systeem af te tappen met de module Harvest (Aftappen) op de touchscreenconsole. **Opmerking:** U moet de module Harvest (Aftappen) configureren voor u deze gebruikt. Zie paragraaf 3.5.8 voor informatie over het configureren van de module Harvest (Aftappen).

 Zoek de module Harvest (Aftappen) op het startscherm van de touchscreenconsole op. Raak de module aan om het scherm Harvest Settings (Aftapinstellingen) te openen en selecteer vervolgens de knop 'Harvest Control' (Aftappen reguleren).

- 2. In het geopende scherm wordt de gebruiker gevraagd een aftapmethode te selecteren: met zwaartekracht of met een pomp.
 - Als u Gravity (Zwaartekracht) selecteert: Voer het vloeistofdoel in. Raak vervolgens 'Done' (Gereed) aan.
 - Als u Pump (Pomp) selecteert: Voer het vloeistofdoel in door de hoeveelheid in te voeren die van de huidige waarde wilt wegnemen. Voer de stroomsnelheid voor fase 1 in, het percentage afgetapt vóór fase 2 en de stroomsnelheid voor fase 2. Raak vervolgens 'Done' (Gereed) aan.
- Als het systeem aangesloten knijpventielen bevat, wordt u in het volgende geopende scherm gevraagd te controleren of de het knijpventiel voor aftappen is gesloten. Als het knijpventiel open is, verplaatst u de schuifknop van 'Open' naar 'Closed' (Gesloten). Raak vervolgens 'Start Harvest' (Aftappen starten) aan.
- 4. Nadat u 'Start Harvest' (Aftappen starten) hebt geselecteerd, worden de pompen ingeschakeld en begint het aftappen van de door de gebruiker gespecificeerde hoeveelheid vloeistof uit het systeem. Het geopende scherm (afbeelding 6.11) toont de huidige fase van het aftapproces de resterende hoeveelheid vloeistof, de huidige massa, de BPC-druk en de vloeistofdruk (vloeistofdruk1, vloeistofdruk2 en delta).
- U kunt te allen tijde onder aan het scherm de knop 'Pause' (Pauzeren) selecteren om het aftapproces tijdelijk te stoppen. Er verschijnt een scherm Harvest Paused (Aftappen gepauzeerd) waarmee u het vullen met vloeistof kunt hervatten of annuleren.



Afbeelding 6.11. Het scherm Harvest Control (Aftappen reguleren).

6.6.2 Handmatig aftappen

Als u geen pompen volgens sectie 3.1 gebruikt, moet u de S.U.M. handmatig aftappen. Volg de onderstaande instructies voor handmatig aftappen.

Handmatig aftappen met een recirculatielus

Als u een recirculatielus gebruikt, bevindt de afvoerleiding zich op toevoegleiding 'Y'. Open de klem op het gedeelte van 30,5 cm (12 inch) van de slang in recirculatielus 'Y' en sluit deze aan op de beoogde overdrachtleiding. Gebruik de peristaltische pomp die al is geïnstalleerd op de recirculatielijn (paragraaf 5.3) om de inhoud van de S.U.M. over te dragen. Zorg dat u de mengmotor stopt voordat u het minimale werkvolume bereikt.

Handmatig aftappen zonder een recirculatielus

Als u geen recirculatielus gebruikt, volgt u de onderstaande stappen om handmatig vloeistof uit de S.U.M. BPC af te tappen.

- 1. Sluit de afvoerslangenset aan de onderkant aan op de beoogde overdrachtslijn. **Opmerking:** Voor steriele aansluitingen moet dit aseptisch worden gedaan.
- 2. Open de klem op de onderste aftappoort.
- 3. Begin de BPC af tappen met behulp van een peristaltische pomp. Zorg dat u de mengmotor stopt voordat de vloeistof in de BPC het minimale werkvolume bereikt.
- Wanneer er ongeveer 3-5 liter in de BPC resteert, tilt u de BPC uit de bovenkant. Houd tijdens het optillen de onderste afvoerleiding bij de vloer om het aftappen van de laatste liter vloeistof te vergemakkelijken.

6.7 Afsluiten en wegwerpen

6.7.1 De S.U.M. afsluiten

 Nadat de S.U.M. is afgetapt, controleert u of de motor van het roermechanisme is uitgeschakeld en schakelt u de voeding naar de buitenste ondersteuningscontainer uit door de hoofdstroom uit te schakelen.

- Verwijder de aandrijfas en berg deze op. Voer hiervoor de stappen voor montage (paragraaf 5.6) in omgekeerde volgorde uit.
 Opmerking: Wees voorzichtig wanneer u de S.U.M. verwijdert onder een laag plafond om beschadiging van het plafond of plafondlampen te voorkomen (zie de vereisten voor plafondhoogte in hoofdstuk 8).
- Als de hardware van de S.U.M. in contact is gekomen met bijtende materialen, dan spoelt u desbetreffende gebieden af met een beetje water en reinigt u ze volgens de standaard procedures. Zie Hoofdstuk 7 – Onderhoud voor meer informatie.
- 4. Breng losse voorwerpen zoals de aandrijfas en gereedschap terug naar de opslaglocaties om accidentele schade te voorkomen.
- 5. Verwijder de BPC van de buitenste ondersteuningscontainer en werp de BPC weg volgens de voorschriften van uw faciliteit. Alle contactmaterialen van het product kunnen in een geschikte afvalbak of verbrandingsinstallatie worden weggedaan. Als de herbruikbare lagerpoort is gebruikt voor open-top mengen, bewaart u deze voor toekomstig gebruik.

6.7.2 De touchscreenconsole buiten bedrijf stellen

- 1. Zet de netstroomschakelaar en de interne wisselstroomschakelaar uit.
- 2. Maak de voedingskabel los.
- 3. Zet de uitvoerschakelaar voor wisselstroom uit.
- 4. Maak de M12-voedingskabel en de motorkabel los. Maak vervolgens alle andere M12-kabels los.
- 5. Maak de optionele externe E-Stop los, indien aanwezig.
- 6. Maak de bevestigingsschroef voorzichtig los en demonteer de eenheid.
- 7. Bewaar de eenheid na verwijdering in een geschikte bergruimte.

6.7.3 Afvoerinformatie

Voer gebruikte BPC's af volgens de voorschriften van uw faciliteit. Alle contactmaterialen van het product kunnen in een geschikte afvalbak of verbrandingsinstallatie worden weggedaan. Producten met het onderstaande symbool voldoen aan richtlijn 2002/96/EC inzake afgedankte elektrische en elektronische apparatuur (AEEA) van de Europese Unie.



Afbeelding 6.12. AEEAsymbool

Als u producten van Thermo Scientific gebruikt, u zich in Europa bevindt en onze producten moet laten recyclen conform de AEEArichtlijn, moet uw product op of na 13 augustus 2005 zijn gekocht en zijn gemarkeerd met het AEEA-symbool. Producten van Thermo Scientific blijven voldoen aan de vereisten van alle relevante richtlijnen, met inbegrip van die van de Richtlijn betreffende de beperking van gevaarlijke stoffen (RoHS) indien van toepassing. Al onze producten waarop de RoHS betrekking heeft of voor het eerst betrekking heeft ten gevolge van de opname in de herziene richtlijn voor productcategorieën 8, 9 of (de nieuwe categorie) 11, wijzigingen in de uitsluitingen of wijzigingen in uitgezonderde toepassingen, zullen vanaf de eerste vervaldatum voldoen aan deze nieuwe vereisten of dit mogelijk maken voor een eindproduct waarin ze worden opgenomen.

Producten die als biologisch gevaarlijk worden beschouwd of medisch besmet zijn, mogen niet via dit programma worden gerecycled. Ze moeten als biologisch gevaarlijk afval worden behandeld en volgens uw lokale voorschriften worden opgeruimd.

Onderhoud en probleemoplossing

Inhoudsopgave

- 7.1 Onderhoud
- 7.2 Probeemoplossing en veelgestelde vragen

7.1 Richtlijnen voor onderhoud

7.1.1 Richtlijnen voor routineonderhoud

Omgevingsomstandigheden, gebruiksparameters en de naleving door de gebruiker van standaard operationele procedures zoals uiteengezet in deze handleiding, kunnen veel invloed hebben op de nuttige levensduur van de S.U.M.-hardware. Er is in het ontwerp van de wegwerpbare constructie van de S.U.M. rekening gehouden met sterk slijtende items die veel voorkomen in conventionele systemen, zoals lagers, afdichtingen, O-ringen en sterilisatiekleppen. Dit heeft een mengsysteem opgeleverd dat inherent robuust is en weinig routineonderhoud vereist. De volgende richtlijnen voor routineonderhoud zijn gebaseerd op standaard gebruiksomstandigheden, zoals gedefinieerd in deze gebruikershandleiding.

Neem tussen het mengen van batches de tijd om de buitenkant van de S.U.M. te reinigen. Dit verbetert het uiterlijk en de algemene levensduur van het hardwaresysteem. Tussen runs kunnen de buitenste ondersteuningscontainer, de aandrijfas, de aandrijving van de menger en de touchscreenconsole worden afgeveegd met schone doekjes. Stalen oppervlakken op de buitenste ondersteuningscontainer kunnen ook gereinigd worden met een reinigingsmiddel voor roestvrij staal. **Opmerking:** Gebruik alleen schone, droge doekjes om het touchscreen van de touchscreenconsole af te vegen.

De S.U.M.-hardware is gemaakt in overeenstemming met de IP-54classificatie voor bescherming tegen binnendringen en kan worden gereinigd volgens de standaardreinigingsprocedures voor laboratoria. Zorg dat alle elektrische aansluitingen zijn losgekoppeld en dat elektriciteitskasten goed gesloten zijn. De unit moet volledig kunnen drogen voordat deze opnieuw in gebruik wordt genomen.

7.1.2 Softwareonderhoud van touchscreenconsole

Wanneer beschikbaar moeten gebruikers software-updates op de touchscreenconsole installeren. U doet dit door de USB-stick met de laatste software-update in de USB-poort van de touchscreenconsole te plaatsen. Selecteer het instellingenpictogram op het startscherm. Klik daarna achtereenvolgens op 'Maintenance' (Onderhoud) en 'Check for updates' (Controleren op updates).

7.1.3 Richtlijnen voor preventief onderhoud

Volg de onderstaande richtlijnen voor preventief onderhoud om betrouwbaar systeemgebruik te garanderen. Aan slijtage onderhevige onderdelen moeten voor en na gebruik op het oog worden geïnspecteerd. Zie de tekeningen in het ETP voor informatie over vervangingsonderdelen. **Opmerking:** De herbruikbare lagerpoort van de S.U.M. moet jaarlijks worden vervangen.

Aandrijfmotor

De aandrijfmotor is een inductiemotor van industriële kwaliteit met een permanent afgedichte en gesmeerde tandwielkast. De aandrijfmotor kan minimaal 10.000 uur met continue belasting worden gebruikt voordat deze moet worden vervangen.

Aandrijfas-eenheid

De aandrijfas slijt bij gebruik enigszins en moet na elke run op het oog worden geïnspecteerd. Coat de uit meerdere segmenten bestaande aandrijfas lichtjes met voor voeding goedgekeurde anti-seize om te helpen bij het maken van de verbindingen van de aandrijfas. De aandrijfas-eenheid moet meestal na één jaar gebruik worden vervangen. Zie voor speciale omstandigheden de slijtagespecificaties van de kop van de aandrijfas in tabel 7.1.

Tabel 7.1 toont de minimale diameters voor de zeskantbus van de aandrijfas. De diameters worden gemeten op de breedste plaats over de punten. Vervang een versleten aandrijfaskop-eenheid wanneer de diameter van de zeskantbus van de aandrijfas op de breedste plek gelijk is aan of kleiner is dan de meting tussen de verschillende punten. Zie afbeelding 7.1 voor de meetlocatie.

Tabel 7.1. Diameters van de zeskantbus van de aandrijfaskop.

Grootte van de S.U.M. Minimale diameter van de zes van de aandrijfas	
50, 100, 200, 500 en 1.000 l	14,4 mm (0,566 inch)
2.000 I	20,8 mm (0,820 inch)



Afbeelding 7.1. Aandrijfaskop-eenheid.

7.2 Problemen met de BPC oplossen

Probleem: Er staat teveel spanning op de film in de buurt van de lagerpoort van de geladen BPC.

Oplossing: Laad zo mogelijk de BPC opnieuw en volg daarbij nauwgezet de BPC-laadinstructies in deze publicatie.

- Zorg voor extra film in het gebied van de BPC dat zich het dichtst bij de lagerpoort bevindt wanneer u de BPC in de tank uitlijnt.
- Vul pas vloeistof toe wanneer de BPC juist is aangebracht in de tank.
- Controleer of de zwarte bumper in het motorsteunblok aanwezig is.

Probleem: Ik ben niet vertrouwd met het gebruik van Pall Kleenpakconnectoren en ben bezorgd over het maken van aseptische verbindingen.

Oplossing: Lees de instructies voor Pall Kleenpak-connectoren in paragraaf 5.7 van deze handleiding voordat u aseptische verbindingen gaat maken.

- Wanneer u een verbinding maakt, moet u de status van de vier externe vergrendelclips beoordelen en controleren of ze goed dicht zitten. U moet bij alle vier de clips een duidelijke klik horen wanneer u de connectoren dicht drukt. Zorg altijd dat de vier vergrendelclips volledig verbonden zijn voor de mannelijke/ vrouwelijke aansluiting voordat u de papieren stroken verwijdert.
- Een veelvoorkomende oorzaak van lekkende Kleenpakconnectoren is een fout in de laatste stap waarbij de tapse cilinders van de mannelijke/vrouwelijke connector goed moeten worden aangebracht. In de mannelijke connector bevindt zich een serie concentrische ringen (0,8 mm/0,3 inch voor de zwarte O-ring). Controleer op het oog of de vier interne clips zich op de

laatste set ringen bevinden. Plaats de connectorflenzen tussen de wijsvinger en duim van uw beide handen en knijp totdat alles goed vastzit.

Probleem: Ik heb voorafgaand aan het vullen met vloeistof geen pH-sonde aangebracht. Kan ik onder die omstandigheden toch nog een steriele aansluiting tot stand brengen?

Oplossing: U kunt nog steeds een steriele aansluiting tot stand brengen nadat u de BPC hebt gevuld. Dit is op voorwaarde dat de klemmen op de sondepoorten van de Kleenpak-connector voorafgaand aan het vullen met vloeistof waren gesloten. De Kleenpak-connectoren moeten droog zijn als u de sondeeenheden wilt aansluiten. Wanneer er al media in de S.U.M. aanwezig is, volgt u de procedures om een sonde in te brengen in paragraaf 5.8 van deze handleiding. Er kan wat vloeistof in de balgen binnendringen wanneer de sonde wordt ingebracht in een BPC die al met vloeistof is gevuld.

7.3 Problemen met de touchscreenconsole oplossen

Zie tabel 7.2 om problemen met de touchscreenconsole op te lossen en potentiële oorzaken te achterhalen. Voor verdere hulp om technische problemen met de touchscreenconsole op te lossen kunt u bellen naar +1 435 792 8500 (Verenigde Staten) of +44 (1) 670 734 093 (Europa, V.K.), of contact opnemen met: customerservice.bioprocessing@thermofisher.com



VOORZICHTIG: Open de touchscreenconsole niet en probeer deze niet zelf te repareren. De eenheid mag alleen worden gerepareerd door servicepersoneel van Thermo Scientific. Reparaties die niet door medewerkers van Thermo Fisher Scientific zijn uitgevoerd, maken de systeemgarantie ongeldig.

Tabel 7.2. Problemen met de touchscreenconso	le oplossen ei	n potentiële oorzaken achterhalen.
--	----------------	------------------------------------

Probleem	Potentiële oorza(a)k(en)
De touchscreenconsole kan niet worden ingeschakeld	De stekker of het stopcontact is defectEen zekering (stroomonderbreker) is doorgeslagen
De touchscreenconsole loopt van op het opstartscherm	 Als dit periodiek gebeurt, kan dit te wijten zijn aan een leesfout, waarvoor opnieuw opstarten vereist is. De firmware is beschadigd of er is een hardwarestoring
De motor start niet	Probleem met de communicatiekabel van de motor
Het toerental (RPM) van de motor valt buiten de tolerantiegrenzen	Dit probleem doet zich niet voor bij gelijkstroommotoren: een optische sensor zorgt dat het toerental binnen de tolerantiegrenzen blijft.
De massa-aflezing wordt niet weergegeven	Probleem met laadcel/laadcelkabelProbleem met optelblok voor laadcellen
De massa-aflezingen zijn onnauwkeurig	De vatgrootte is onjuistDe laadcelkalibratie is onnauwkeurig
De pH-aflezingen worden niet weergegeven of de pH- zender herkent de sensor niet.	Probleem met pH-verbindingProbleem met pH-kaart in de touchscreenconsole
De pH-aflezingen zijn onnauwkeurig	 pH-aflezingen worden altijd weergegeven, zelfs zonder aangesloten sensor of kabel Probleem met pH-kalibratie
De geleidbaarheidsaflezingen worden niet weergegeven	Probleem met de geleidbaarheidssensorProbleem met de geleidbaarheidsverbinding
De geleidbaarheidsaflezingen zijn onnauwkeurig	Probleem met de geleidbaarheidskalibratieProbleem met de geleidbaarheidssensor
De BPC-druk wordt niet afgelezen	 De druksensor in de BPC werkt niet Verbindingsprobleem tussen een druksensor en de touchscreenconsole
De aflezingen van de BPC-druk zijn onnauwkeurig	 Probleem met de kalibratie van de druksensor Verbindingsprobleem tussen een druksensor en de touchscreenconsole
De BPC-drukmodule pompt geen lucht in de BPC nadat op 'Start' (Starten) is gedrukt	Probleem met luchttoevoer in de touchscreenconsoleProbleem met de pneumatische magneetkleppen

Tabel 7.3. Problemen met de touchscreenconsole oplossen en potentiële oorzaken achterhalen (vervolg).

Probleem	Potentiële oorza(a)k(en)
De touchscreenconsole regelt de TCU niet	Probleem met TCU-parametersProbleem met TCU-kabelsets
Een onderbreker van de touchscreenconsole slaat binnen een periode van vier uur meerdere keren door	 Inadequate stroomvoorziening: zorg dat er geen andere apparaten op hetzelfde stopcontact zijn aangesloten. Er wordt teveel stroom verbruikt door (al dan niet defecte) apparaten die op hulpstopcontacten zijn aangesloten. Verwijder het of de apparaten om de defecte component te vinden
Het HMI-scherm van de touchscreenconsole reageert niet	Probleem met de softwareProbleem met de elektrische hardware
De touchscreenconsole herkent een USB-stick niet	De USB-stick heeft geen FAT-indelingProbleem met elektronische hardware
De touchscreenconsole regelt een knijpventiel niet	 Onjuiste feedback- en controleaansluiting van knijpventiel Probleem met persluchtingang
De touchscreenconsole regelt een extern apparaat niet (uitgang van 4-20 mA)	 Probleem met parameters op externe component Probleem met kabelset
De touchscreenconsole herkent een hulpingang niet (ingang van 4-20 mA)	 Probleem met externe toevoer van 4-20 mA Probleem met kabelset
Een hulponderbreker slaat binnen een periode van vier uur meerdere keren door	Overbelasting van hulpstekkersProbleem met onderbreker
De temperatuuraflezingen worden niet weergegeven of herkennen een sensor niet	Probleem met RTDProbleem met kabelset
De temperatuuraflezingen zijn onnauwkeurig	Probleem met RTDDe temperatuurkalibratie is onnauwkeurig
Het digitale communicatieprotocol werkt niet	 Probleem met de software Probleem met kabelset
De externe E-Stop wordt niet herkend	Probleem met kabelsetProbleem met het circuit van de E-Stop
De zoemer van de E-Stop kan niet worden uitgeschakeld	Probleem met het resetcircuit van de E-StopProbleem met zoemer
De zoemer van de E-Stop gaat niet aan wanneer er op de knop van de E-Stop wordt gedrukt	Probleem met zoemerProbleem met het circuit van de E-Stop
De hulpcontacten werken niet	 Een hulponderbreker is mogelijk doorgeslagen Probleem met netsnoer
Waarschuwing voor oververhitting	Probleem met uitlaatventilatorProbleem met inlaatventilator
HMI reageert langzaam of loopt achter	Probleem met firmware
De gelijkstroommotor is oververhit	Probleem met koelventilatorProbleem met bedrading

0

Specificaties en informatie over onderdelen

Inhoudsopgave

- 8.1 Hardware-kenmerken
- 8.2 Hardwarespecificaties
- 8.3 Specificaties van de touchscreenconsole
- 8.4 BPC-specificaties
- 8.5 Accessoires en opties

8.1 Hardware-kenmerken

8.1.1 Ontwerpkenmerken van S.U.M.'s van 50, 100 en 200 l



Afbeelding 8.1. S.U.M. met mantel, 200 I (vooraanzicht).

- 1. Powdertainer-arm (optioneel)
- 2. Mengsysteem met afscherming
- 3. Steun van lagerpoort met klem
- 4. Mixermotor
- 5. Aandrijfas, opgeborgen
- 6. Planken (optioneel)
- 7. Kijkvenster voor vloeistof
- 8. Touchscreenconsole
- 9. Sondecliphanger
- 10. Vensters voor sondetoegang
- 11. Wagen
- 12. Kabelbeheerhaken

Afbeelding 8.2. S.U.M. met mantel, 200 I (achteraanzicht).

- Standaard gereedschapsset: Vierkante momentsleutel van 10 mm (3/8 inch) x 16,9 Nm (150 in-lb.) en vergrendelsleutel voor laadcellen en motorkap
- 14. Gegolfde mantel 0,95 cm (3/8 inch)
- 15. Buitenste ondersteuningscontainer van roestvrij staal
- 16. Handgrepen met clips voor kabelbeheer
- 17. Mand (optioneel)
- 18. Openingen/pennen onderaan voor bevestiging en uitlijning BPC
- 19. Aftapklep (alleen op modellen met mantel)
- 20. 3,81 cm (1,5 inch) poorten voor aansluiting met drie klemmen voor watertoevoer/-afvoer (alleen op modellen met mantel)
- 21. Wielen (2 draaiend, 2 vast)

8.1.2 Ontwerpkenmerken van S.U.M.'s van 500 en 1.000 l



Afbeelding 8.3. S.U.M. met mantel, 500 I (vooraanzicht).

- 1. Powdertainer-arm (optioneel)
- 2. Mengeenheid met veiligheidsdeksel
- 3. Steun van lagerpoort met klem
- 4. Mixermotor
- 5. Aandrijfas, opgeborgen
- 6. Planken (optioneel)
- 7. Deur om BPC te laden, met kijkvensters voor vloeistof
- 8. Touchscreenconsole
- 9. Sondecliphanger
- 10. Vensters voor sondetoegang
- 11. Wagen
- 12. Kabelbeheerhaken
- Standaard gereedschapsset: Vierkante momentsleutel van 10 mm (3/8 inch) x 16,9 Nm (150 in-lb.) en vergrendelsleutel voor laadcellen en motorkap



Afbeelding 8.4. S.U.M. met mantel, 500 I (achteraanzicht).

- 14. Gegolfde mantel 0,95 cm (3/8 inch)
- 15. Buitenste ondersteuningscontainer van roestvrij staal
- 16. Handgrepen met clips voor kabelbeheer
- 17. Mand (optioneel)
- 18. Uitsparing met verwijderbare beugel om BPC te laden
- 19. Openingen/pennen onderaan voor bevestiging en uitlijning BPC
- 20. Aftapklep (alleen op modellen met mantel)
- 21. 3,81 cm (1,5 inch) poorten voor aansluiting met drie klemmen voor watertoevoer/-afvoer (alleen op modellen met mantel)
- 22. Wielen (2 draaiend, 2 vast)

8.1.3 Ontwerpkenmerken van S.U.M.'s van 2.000 l



Afbeelding 8.5. S.U.M. met mantel, 2.000 I (vooraanzicht).

- 1. Powdertainer-arm (optioneel)
- 2. Mengeenheid met veiligheidsdeksel
- 3. Steun van lagerpoort met klem
- 4. Mixermotor
- 5. Touchscreenconsole
- 6. Aandrijfas, opgeborgen
- 7. Kabelbeheerhaken
- 8. Venster sondetoegang
- 9. Sondecliphanger
- 10. BPC-laaddeur en kijkvensters voor vloeistof
- 11. Mand (optioneel)



Afbeelding 8.6. S.U.M. met mantel, 2.000 I (achteraanzicht).

- 12. Planken (optioneel)
- Standaard gereedschapsset: Vierkante momentsleutel van 10 mm (3/8 inch) x 16,9 Nm (150 in-lb.) en vergrendelsleutel voor laadcellen en motorkap
- 14. Gegolfde mantel 0,95 cm (3/8 inch)
- 15. Buitenste ondersteuningscontainer van roestvrij staal
- 16. Achterdeur (om BPC te laden), met kijkvenster
- 17. Openingen/pennen onderaan voor bevestiging en uitlijning BPC
- 18. Aftapklep (alleen op modellen met mantel)
- 3,81 cm (1,5 inch) poorten voor aansluiting met drie klemmen voor watertoevoer/-afvoeren (alleen op modellen met mantel)
- 20. Stelpoten (3)

8.2 Hardware-specificaties

De volgende tabellen en afbeeldingen bevatten specificatiegegevens voor S.U.M.-systemen van 50, 100, 200, 500, 1.000 en 2.000 l

Opmerking: Afbeeldingen 8.7 tot en met 8.18 tonen mengers met watermantels. Modellen zonder watermantel kunnen iets andere afmetingen hebben. Zie de bij uw eenheid geleverde tekeningen voor de exacte maten.

Tabel 8.1. Specificaties voor S.U.M.'s van 50 I

		Alleen gelijkstroommotoren	
		Mantel	Zonder mantel
	Nominaal werkvolume van vloeistof	50	
	Minimaal werkvolume van vloeistof	10	
	Totaal kamervolume (vloeistof en gas)	80 I	
etrie	Diameter BPC-kamer	34,9 cm (13,75 inch)	
eome	Schouderhoogte BPC-kamer	84,8 cm (33	,4 inch)
gerg	Vloeistofhoogte en nominaal werkvolume	52,1 cm (20	,5 inch)
Men	Vloeistofgeometrie bij verhouding (hoogte/diameter) werkvolume	1,5:1	
	Ophopingsvolume	< 50 r	nl
	Algemene mengergeometrie (verhouding hoogte/diameter)	1,9:1	
	Tankschotten	Geen	
ef	Mengschroef (hoeveelheid x aantal schoepen)	1 x 3	
chro	Schaal van mengschroef (diameter mengschroef/diameter tank)	2/5	
engs	Pitch van mengschroefschoepen (hoek)	45°	
Σ	Mengschroefdiameter	14,6 cm (5,7	75 inch)
	Bereik mengsnelheid	20–356 I	RPM
	Snelheid van tip	15 cm/s (28,5 ft/min) – 27	2 cm/s (535,4 ft/min)
	Mengstroomrichting linksom	Omlaag pompen	
en	Opgeloste hoek roeras	12,5°)
hudd	Verschuiving van middellijn roeras	1,9 cm (0,75 inch)	
Sc	Totale lengte aandrijfas	91,7 cm (36	,1 inch)
	Diameter aandrijfas	1,27 cm (0.	5 inch)
	Buitendiameter plastic omhulsel aandrijfas	2,54 cm (1	inch)
	Vrije ruimte mengschroef vanaf tankbodem	11,75 cm (4,63 inch)	

Tabel 8.2. Specificaties voor S.U.M.'s van 50 I (vervolg).

		Alleen gelijkstroommotoren	
		Mantel	Zonder mantel
for	Motoraandrijving roermechanisme (type, spanning, fase)	Borstelloos, 48 VDC	
	Nominaal motorvermogen	400 W (0,536 pk) Motor: 48 V, 6,5 A	
Moi	Nominaal motorkoppel	8,6 Nm (76	in-lb.)
	Tandwielreductie	7,5:1	
	Methodes voor motorcommunicatie	Op de touchscreenconsole via communicatiepoorten	
antel	Manteloppervlak: vol/half volume	0,38 m ² (4,1 ft ²) / 0,32 m ² (3,4 ft ²)	-
stofn	Mantelvolume	2 I (0,53 gal.)	-
loeis	Manteldebiet bij 3,4 bar (50 psi)	99 l/min (26,4 gal/min)	-
>	Procesverbinding	Hygiënische drieklem 1 inch	-
ė	TCU-model: maximale verwarming/koeling	TF2500: 2800/2500 W	-
tuurr ing	Geschatte opwarmtijd voor vloeistof (5–37°C)	1,2 uur	-
erat uleri	Geschatte afkoeltijd voor vloeistof (37-5°C)	2,7 uur	-
Temp	RTD of thermokoppel, 3,18 mm (1/8 inch) OD	Pt-100 (standaard)	
ainer and)	Totale breedte	87 cm (34,1 inch)	
scont: en/ma	Totale lengte	86 cm (33,9 inch)	
uning: Jank	Algehele hoogte (zonder Powdertainer-arm)	152 cm (59,8 inch)	
erstei nder j	Gewicht droge skid (massa)	164,1 kg (360,8 lb.)	145,1 kg (319,8 lb.)
Ondi (zo	Gewicht natte skid–nominaal werkvolume (massa)	214 kg (470,8 lb.)	195,1 kg (429,8 lb.)
	Vereiste plafondhoogte voor laden van aandrijfas	228,6 cm (90 inch)	
	Nominale voeding	100-120 VAC, 50/60 Hz, enkel, 15 A 220-240 VAC, 50/60 Hz, enkel, 10,4 A	
emeen	pH- en DO-sonde —autoclaveerbaar type (Applisens, Broadley James, Mettler-Toledo)	12 mm diameter x 215–235 mm invoerlengte x 13,5 PG (pijp) schroefdraad	
Alg	Ruisniveau	< 70 dB op 1,5 m	
	Opslagtemperatuur	-25–65°C	
	Relatieve vochtigheid	20-80% zonder (condensatie
S	Bereik van bedrijfstemperatuur	S.U.M.: 2–40 +/- 0,1°C (36–104 +/- 0,2 °F) Gelijkstroommotor 0–40°C	
en nete	Motorsnelheid	20–356 F	RPM
evole aran	Volumebereik	10–50	1
Aanbi uiksp	Maximale statische BPC-druk	0,03 bar (0	,5 psi)
gebr	Maximale BPC-druk tijdens gebruik	0,007 bar (0),1 psi)
0)	Continue bedrijfsduur	21 dagen mengen alleen bij nominaal volume	



Afbeelding 8.7. Afmetingen van S.U.M. van 50 I (vooraanzicht).



Afbeelding 8.8. Afmetingen van S.U.M. van 50 I (bovenaanzicht).

Tabel 8.3. Bestelgegevens voor S.U.M.'s van 50 l.

Beschrijving	Cat.nr.
50 I zonder mantel, gelijkstroommotor, touchscreenconsole, met laadcellen	SUM0050.9001
50 I met mantel, gelijkstroommotor, touchscreenconsole, met laadcellen	SUM0050.9002

Tabel 8.4. Specificaties voor S.U.M.'s van 100 I

		Alleen gelijkstroommotoren	
		Mantel	Zonder mantel
	Nominaal werkvolume van vloeistof	100	
	Minimaal werkvolume van vloeistof	20 I (met complete mengschroefbedekking)	
	Totaal kamervolume (vloeistof en gas)	150 l	
etrie	Diameter BPC-kamer	43,8 cm (17,2	5 inch)
eom	Schouderhoogte BPC-kamer	100,8 cm (39,7 inch)	
gerg	Vloeistofhoogte en nominaal werkvolume	66 cm (26 inch)	
Men	Vloeistofgeometrie bij verhouding (hoogte/diameter) werkvolume	1,5:1	
	Ophopingsvolume	< 50 m	1
	Algemene mengergeometrie (verhouding hoogte/diameter)	1,9:1	
	Tankschotten	Geen	
ef	Mengschroef (hoeveelheid x aantal schoepen)	1 x 3	
chro	Schaal van mengschroef (diameter mengschroef/diameter tank)	1/3	
sbue	Pitch van mengschroefschoepen (hoek)	45°	
Ň	Mengschroefdiameter	14,6 cm (5,75 inch)	
	Bereik mengsnelheid	20–356 RPM	
	Snelheid van tip	22,9 cm/s (45,2 ft/min) – 267,7 cm/s (526,9 ft/min)	
	Mengstroomrichting linksom	Omlaag pompen	
den	Opgeloste hoek roeras	15°	
pnuc	Verschuiving van middellijn roeras	5,08 cm (2	inch)
Š	Totale lengte aandrijfas	104,4 cm (41,1 inch)	
	Diameter aandrijfas	1,27 cm (0.5	inch)
	Buitendiameter plastic omhulsel aandrijfas	2,54 cm (1 inch)	
	Vrije ruimte mengschroef vanaf tankbodem	4,9 cm (1,93	inch)
	Motoraandrijving roermechanisme (type, spanning, fase)	Borstelloos, 4	8 VDC
z	Nominaal motorvermogen	400 W (0,536 pk) Motor: 48 V, 6,5 A	
Mot	Nominaal motorkoppel	8,6 Nm (76 i	n-lb.)
	Tandwielreductie	7,5:1	
	Methodes voor motorcommunicatie	Op de touchscreenconsole via communicatiepoorten	

Tabel 8.5. Specificaties voor S.U.M.'s van 100 I (vervolg).

		Alleen gelijkstroommotoren	
		Mantel	Zonder mantel
ntel	Manteloppervlak: vol/half volume	0,69 m ² (7,4 ft ²) / 0,41 m ² (4,4 ft ²)	-
ofma	Mantelvolume	4 I (1,1 gal.)	-
eist(Manteldebiet bij 3,4 bar (50 psi)	100 l/min (26,4 gal/min)	-
VIO	Procesverbinding	Hygiënische drieklem 1 inch	-
regu-	TCU-model: maximale verwarming/koeling	TF2500: 2800/2500 W	-
ıtuur ring	Geschatte opwarmtijd voor vloeistof (5–37°C)	2 uur	-
ipera lei	Geschatte afkoeltijd voor vloeistof (37–5°C)	5,1 uur	-
Tem	RTD of thermokoppel, 3,18 mm (1/8 inch) OD	Pt-100 (s	tandaard)
ainer ind)	Totale breedte	87 cm (3	4,1 inch)
scont en/m	Totale lengte	92 cm (36,4 inch)	
ining: Jank	Algehele hoogte (zonder Powdertainer-arm)	153 cm (60,2 inch)	
ersteu 1der p	Gewicht droge skid (massa)	172 kg (379,8 lb.)	194,1 kg (426,8 lb.)
0nde (zoi	Gewicht natte skid–nominaal werkvolume (massa)	272,1 kg (600,8 lb.)	294,1 kg (647,8 lb.)
	Vereiste plafondhoogte voor laden van aandrijfas	236,2 cm (93 inch)	
	Nominale voeding	100–120 VAC, 50/60 Hz, enkel, 15 A 220–240 VAC, 50/60 Hz, enkel, 10,4 A	
Jemeen	pH- en DO-sonde —autoclaveerbaar type (Applisens, Broadley James, Mettler-Toledo)	12 mm diameter x 215–235 mm invoerlengte x 13,5 PG (pijp) schroefdraad	
Alĝ	Ruisniveau	< 70 dB	op 1,5 m
	Opslagtemperatuur	-25–65°C	
	Relatieve vochtigheid	20–80% zonde	er condensatie
S	Bereik van bedrijfstemperatuur	S.U.M.: $2-40 \pm 0,1^{\circ}C (36-104 \pm 0,2^{\circ}F)$ Gelijkstroommotor $0-40^{\circ}C$	
en neter	Motorsnelheid	20–35	6 RPM
ievole parar	Volumebereik	20–100 I	
Aanb ruiks	Maximale statische BPC-druk	0,03 bar	(0,5 psi)
gebr	Maximale BPC-druk tijdens gebruik	0,007 ba	r (0,1 psi)
	Continue bedrijfsduur	21 dagen mengen alleen bij nominaal volume	



Afbeelding 8.9. Afmetingen van S.U.M. van 100 I (vooraanzicht).



Afbeelding 8.10. Afmetingen van S.U.M. van 100 I (bovenaanzicht).

Tabel 8.6. Bestelgegevens voor S.U.M.'s van 100 l.

Beschrijving	Cat.nr.
100 I zonder mantel, gelijkstroommotor, touchscreenconsole, met laadcellen	SUM0100.9001
100 I met mantel, gelijkstroommotor, touchscreenconsole, met laadcellen	SUM0100.9002

Tabel 8.7. Specificaties voor S.U.M.'s van 200 I

		Alleen gelijkstroommotoren	
		Mantel	Zonder mantel
	Nominaal werkvolume van vloeistof	200	
	Minimaal werkvolume van vloeistof	40 I (met complete mengschroefbedekking)	
	Totaal kamervolume (vloeistof en gas)	250	
trie	Diameter BPC-kamer	59,7 cm (20 inch)	
ome	Schouderhoogte BPC-kamer	123,4 cm (48,6 inch)	
erge	Vloeistofhoogte en nominaal werkvolume	99 cm (3	39 inch)
Meng	Vloeistofgeometrie bij verhouding (hoogte/diameter) werkvolume	1,5	:1
	Ophopingsvolume	< 50) ml
	Algemene mengergeometrie (verhouding hoogte/diameter)	1,95	5:1
	Tankschotten	Geo	en
-	Mengschroef (hoeveelheid x aantal schoepen)	1 x	3
schroel	Schaal van mengschroef (diameter mengschroef/diameter tank)	2/5	
leng	Pitch van mengschroefschoepen (hoek)	45°	
2	Mengschroefdiameter	20 cm (7,87 inch)	
	Bereik mengsnelheid	20-356 RPM	
	Snelheid van tip	31,4–366,6 cm/s (61,9–721,6 ft/min)	
	Mengstroomrichting linksom	Omlaag pompen	
en	Opgeloste hoek roeras	12,5°	
ppnu	Verschuiving van middellijn roeras	6,4 cm (2	2,5 inch)
Scl	Totale lengte aandrijfas	129,5 cm	(51 inch)
	Diameter aandrijfas	1,27 cm (0.5 inch)
	Buitendiameter plastic omhulsel aandrijfas	2,54 cm	(1 inch)
	Vrije ruimte mengschroef vanaf tankbodem	7,9 cm (3	3,1 inch)
	Motoraandrijving roermechanisme (type, spanning, fase)	Borstelloos	s, 48 VDC
-	Nominaal motorvermogen	400 W (0,536 pk) Motor: 48 V, 6,5 A	
Mot	Nominaal motorkoppel	8,6 Nm (7	76 in-lb.)
-	Tandwielreductie	7,5:1	
	Methodes voor motorcommunicatie	Op de touchscreenconsole via communicatiepoorten	

Tabel 8.8. Specificaties voor S.U.M.'s van 200 I (vervolg).

		Alleen gelijkstroommotoren	
		Mantel	Zonder mantel
tel	Manteloppervlak: vol/half volume	1,2 m² (13 ft²) / 0,69 m² (7,4 ft²)	-
iman	Mantelvolume	6,5 I (1,7 gal.)	-
oeistof	Manteldebiet bij 3,4 bar (50 psi)	99,4 l/min (26,3 gal/min)	-
N	Procesverbinding	Hygiënische drieklem 1 inch	-
rre-	TCU-model: maximale verwarming/koeling	TF2500: 2800/2500 W	-
atuur sring	Geschatte opwarmtijd voor vloeistof (5-37°C)	3,4 uur	-
gule	Geschatte afkoeltijd voor vloeistof (37–5°C)	6,8 uur	-
Ten	RTD of thermokoppel, 3,18 mm (1/8 inch) OD	Pt-100 (s	tandaard)
ai-	Totale breedte	90 cm (3	5,4 inch)
scont anke	Totale lengte	107 cm (42 inch)	
unings der pl nand)	Algehele hoogte (zonder Powdertainer-arm)	174 cm (68,5 inch)	
lerstei r (zon n	Gewicht droge skid (massa)	259,1 kg (570,8 lb.)	230,1 kg (507,8 lb.)
Ond	Gewicht natte skid–nominaal werkvolume (massa)	459,1 kg (1011,8 lb.)	430,1 kg (948,8 lb.)
	Vereiste plafondhoogte voor laden van standaard aandrijfas	281,9 cm (111 inch)	
	Vereiste plafondhoogte voor laden van optionele 2-delige aandrijfas	242,3 cm (95,4 inch)	
een	Nominale voeding	100–120 VAC, 50/60 Hz, enkel, 15 A 220–240 VAC, 50/60 Hz, enkel, 10,4 A	
gem	Opslagtemperatuur	-25–	65°C
AI	Relatieve vochtigheid	20–80% zonder condensatie	
	pH- en DO-sonde —autoclaveerbaar type (Applisens, Broadley James, Mettler-Toledo)	12 mm diameter x 215–235 mm invoerlengte x 13,5 PG (pijp) schroefdraad	
	Ruisniveau	< 70 dB op 1,5 m	
irs	Bereik van bedrijfstemperatuur	S.U.M.: $2-40 \pm 0,1^{\circ}C$ ($36-104 \pm 0,2^{\circ}F$) Gelijkstroommotor $0-40^{\circ}C$	
en mete	Motorsnelheid	20–356 RPM	
evol	Volumebereik	40–200 I	
Aanb uiksp	Maximale statische BPC-druk	0,03 bar	(0,5 psi)
gebri	Maximale BPC-druk tijdens gebruik	0,007 ba	r (0,1 psi)
	Continue bedrijfsduur	21 dagen mengen alleen bij nominaal volume	



Afbeelding 8.11. Afmetingen van S.U.M. van 200 I (vooraanzicht).



Afbeelding 8.12. Afmetingen van S.U.M. van 200 l (bovenaanzicht).

Tabel 8.9. Bestelgegevens voor S.U.M.'s van 200 l.

Beschrijving	Cat.nr.	
200 I zonder mantel, gelijkstroommotor, touchscreenconsole, met laadcellen	SUM0200.9001	
200 I met mantel, gelijkstroommotor, touchscreenconsole, met laadcellen	SUM0200.9002	

Tabel 8.10. Specificaties voor S.U.M.'s van 500 I

		Alleen gelijkstroommotoren	
		Mantel	Zonder mantel
eometrie	Nominaal werkvolume van vloeistof	500 I	
	Minimaal werkvolume van vloeistof	100 I (met complete mengschroefbedekking)	
	Totaal kamervolume (vloeistof en gas)	660	
	Diameter BPC-kamer	75,56 cm (29,75 inch)	
	Schouderhoogte BPC-kamer	146 cm (57,5 inch)	
gerg	Vloeistofhoogte en nominaal werkvolume	111,8 cm (44 inch)	
Menç	Vloeistofgeometrie bij verhouding (hoogte/diameter) werkvolume	1,5:	1
	Ophopingsvolume	< 100) ml
	Algemene mengergeometrie (verhouding hoogte/diameter)	1,7:1	
	Tankschotten	Geen	
ef	Mengschroef (hoeveelheid x aantal schoepen)	1 x 3	
chro	Schaal van mengschroef (diameter mengschroef/diameter tank)	9/3	4
Mengs	Pitch van mengschroefschoepen (hoek)	45°	
	Mengschroefdiameter	20 cm (7,87 inch)	
	Bereik mengsnelheid	20–356 RPM	
	Snelheid van tip	31,4–366,6 cm/s (61,9–721,6 ft/min)	
	Mengstroomrichting linksom	Omlaag pompen	
len	Opgeloste hoek roeras	20°	
pnqq	Verschuiving van middellijn roeras	10,56 cm (4	l,16 inch)
Sc	Totale lengte aandrijfas	152,4 cm ((60 inch)
	Diameter aandrijfas	1,27 cm (0.5 inch)	
	Buitendiameter plastic omhulsel aandrijfas	2,54 cm (1 inch)	
	Vrije ruimte mengschroef vanaf tankbodem	7,52 cm (2,96 inch)	
Motor	Motoraandrijving roermechanisme (type, spanning, fase)	Borstelloos, 48 VDC	
	Nominaal motorvermogen	400 W (0, Motor: 48	536 pk) V, 6,5 A
	Nominaal motorkoppel	-	
	Tandwielreductie	7,5:	1
	Methodes voor motorcommunicatie	Op de touchscreenconsole via communicatiepoorten	

Tabel 8.11. Specificaties voor S.U.M.'s van 500 I (vervolg).

		Alleen gelijkstroommotoren	
		Mantel	Zonder mantel
Vloeistofmantel	Manteloppervlak: vol/half volume	2,2 m² (23,9 ft²) / 1,4 m² (14,8 ft²)	-
	Mantelvolume	11 I (2,9 gal.)	-
	Manteldebiet bij 3,4 bar (50 psi)	96,8 l/min (25,5 gal/min)	-
	Procesverbinding	Hygiënische drieklem 1 inch	-
-e-	TCU-model: maximale verwarming/koeling	TF10000: 6100/10000 W	-
atuur ring	Geschatte opwarmtijd voor vloeistof (5–37°C)	2,6 uur	-
ipera gule	Geschatte afkoeltijd voor vloeistof (37–5°C)	3,7 uur	-
Tem	RTD of thermokoppel, 3,18 mm (1/8 inch) OD	Pt-100 (standaard)	
container n/mand)	Totale breedte	118 cm (46,	2 inch)
	Totale lengte	132 cm (51,8 inch)	
lnings Ianke	Algehele hoogte (zonder Powdertainer-arm)	197 cm (77,7 inch)	
ersteu nder p	Gewicht droge skid (massa)	424,1 kg (935,8 lb.)	326,1 kg (718,8 lb.)
Onde (zor	Gewicht natte skid–nominaal werkvolume (massa)	924,1 kg (2037,8 lb.)	826,1 kg (1820,8 lb.)
	Vereiste plafondhoogte voor laden van standaard aandrijfas	320 cm (126 inch)	
	Vereiste plafondhoogte voor laden van optionele 2-delige aandrijfas	275,6 cm (108,5 inch)	
een	Nominale stroomvoorziening	100–120 VAC, 50/60 Hz, enkel, 15 A 220–240 VAC, 50/60 Hz, enkel, 10,4 A	
Algem	pH- en DO-sonde —autoclaveerbaar type (Applisens, Broadley James, Mettler-Toledo)	12 mm diameter x 215–235 mm invoerlengte x 13,5 PG (pijp) schroefdraad	
	Ruisniveau	< 70 dB op 1,5 m	
	Opslagtemperatuur	-25–65°C	
	Relatieve vochtigheid	20–80% zonder condensatie	
ſS	Bereik van bedrijfstemperatuur	S.U.M.: $2-40 \pm 0,1^{\circ}C$ (36–104 ± 0,2 °F) Gelijkstroommotor 0–40°C	
en nete	Motorsnelheid	20–356 RPM	
anbevole iiksparan	Volumebereik	100–500 I	
	Maximale statische BPC-druk	0,03 bar (0,5 psi)	
jebru	Maximale BPC-druk tijdens gebruik	0,007 bar (0,1 psi)	
6)	Continue bedrijfsduur	21 dagen mengen alleen	bij nominaal volume



Afbeelding 8.13. Afmetingen van S.U.M. van 500 I (vooraanzicht).



Afbeelding 8.14. Afmetingen van S.U.M. van 500 I (bovenaanzicht).

Tabel 8.12. Bestelgegevens voor S.U.M.'s van 500 I.

Beschrijving	Cat.nr.
500 I zonder mantel, gelijkstroommotor, touchscreenconsole, met laadcellen	SUM0500.9001
500 I met mantel, gelijkstroommotor, touchscreenconsole, met laadcellen	SUM0500.9002

Tabel 8.13. Specificaties voor S.U.M.'s van 1.000 I

		Alleen gelijkstroommotoren	
		Mantel	Zonder mantel
Mengergeometrie	Nominaal werkvolume van vloeistof	1.000 l	
	Minimaal werkvolume van vloeistof	200 I (met complete mengschroefbedekking)	
	Totaal kamervolume (vloeistof en gas)	1.380 I	
	Diameter BPC-kamer	105,4 cm (41,52 inch)	
	Schouderhoogte BPC-kamer	157,2 cm (61,9 inch)	
	Vloeistofhoogte en nominaal werkvolume	115,6 cm (4	45,5 inch)
	Vloeistofgeometrie bij verhouding (hoogte/diameter) werkvolume	1,1:	:1
	Ophopingsvolume	< 300) ml
	Algemene mengergeometrie (verhouding hoogte/diameter)	1,2:1	
	Tankschotten	Geen	
ef	Mengschroef (hoeveelheid x aantal schoepen)	1 x 3	
Mengschroe	Schaal van mengschroef (diameter mengschroef/diameter tank)	8/2	5
	Pitch van mengschroefschoepen (hoek)	45°	
	Mengschroefdiameter	20 cm (7,87 inch)	
	Bereik mengsnelheid	20-356 RPM	
	Snelheid van tip	31,4-366,6 cm/s (61,9-721,6 ft/min)	
	Mengstroomrichting linksom	Omlaag pompen	
len	Opgeloste hoek roeras	22	0
ppnu	Verschuiving van middellijn roeras	12,7 cm (5 inch)	
Sc	Totale lengte aandrijfas	152,4 cm (60 inch)	
	Diameter aandrijfas	1,27 cm (0.5 inch)	
	Buitendiameter plastic omhulsel aandrijfas	2,54 cm (1 inch)	
	Vrije ruimte mengschroef vanaf tankbodem	11,81 cm (4,65 inch)	
Motor	Motoraandrijving roermechanisme (type, spanning, fase)	Borstelloos	, 48 VDC
	Nominaal motorvermogen	400 W (0, Motor: 48	536 pk) V, 6,5 A
	Nominaal motorkoppel	8,6 Nm (7	6 in-lb.)
	Tandwielreductie	7,5	:1
	Methodes voor motorcommunicatie	Op de touchscre communica	enconsole via tiepoorten
Tabel 8.14. Specificaties voor S.U.M.'s van 1.000 I (vervolg).

		Alleen gelijkstroommotoren		
		Mantel	Zonder mantel	
ntel	Manteloppervlak: vol/half volume	3,5 m² (37,6 ft²) / 2,1 m² (22,7 ft²)	-	
fmai	Mantelvolume	17,5 I (4,6 gal.)	-	
loeisto/	Manteldebiet bij 3,4 bar (50 psi)	102 l/min (27 gal/min)	-	
>	Procesverbinding	Hygiënische drieklem 1 inch	-	
're-	TCU-model: maximale verwarming/koeling	TF24000: 22500/24000 W	-	
atuur ring	Geschatte opwarmtijd voor vloeistof (5–37°C)	1,6 uur	-	
ipera gule	Geschatte afkoeltijd voor vloeistof (37–5°C)	2,3 uur	-	
Tem	RTD of thermokoppel, 3,18 mm (1/8 inch) OD	(1/8 inch) OD Pt-100 (standaard)		
iner nd)	Totale breedte	144 cm (56,4	l inch)	
conta n/mai	Totale lengte	158 cm (62,3	3 inch)	
nings lanke	Algehele hoogte (zonder Powdertainer-arm)	200 cm (78,7 inch)		
rsteu der p	Gewicht droge skid (massa)	559,1 kg (1232,8 lb.)	439 kg (967,8 lb.)	
Onde (zon	Gewicht natte skid–nominaal werkvolume (massa)	1559,1 kg (3437,8 lb.)	1439,1 kg (3172,8 lb.)	
	Vereiste plafondhoogte voor laden van standaard aandrijfas	320 cm (126 inch)		
	Vereiste plafondhoogte voor laden van optionele 2-delige aandrijfas	277,6 cm (109,3 inch)		
leen	Nominale voeding	100–120 VAC, 50/60 Hz, enkel, 15 A 200–240 VAC, 50/60 Hz, enkel, 10,4 A		
lgem	Ruisniveau	< 70 dB op ⁻	1,5 m	
Α	pH- en DO-sonde —autoclaveerbaar type (Applisens, Broadley James, Mettler-Toledo)	12 mm diameter x 215–23 13,5 PG (pijp) sch	5 mm invoerlengte x nroefdraad	
	Opslagtemperatuur	-25–65°	С	
	Relatieve vochtigheid	20–80% zonder condensatie		
ľS	Bereik van bedrijfstemperatuur	S.U.M.: 2–40 ± 0,1°C (36–104 ± 0,2 °F) Gelijkstroommotor 0–40°C		
en mete	Motorsnelheid	20-356 RPM		
evol	Volumebereik	200–1.000 l		
Aanb uiksp	Maximale statische BPC-druk	0,03 bar (0,5	ō psi)	
) Jebri	Maximale BPC-druk tijdens gebruik	0,007 bar (0,1 psi)		
	Continue bedrijfsduur	21 dagen mengen alleen b	ij nominaal volume	



Afbeelding 8.15. Afmetingen van S.U.M. van 1.000 I (vooraanzicht).



Afbeelding 8.16. Afmetingen van S.U.M. van 1.000 I (bovenaanzicht).

Tabel 8.15. Bestelgegevens voor S.U.M.'s van 1.000 l.

Beschrijving	Cat.nr.
1.000 I zonder mantel, gelijkstroommotor, touchscreenconsole, met laadcellen	SUM1000.9001
1.000 I met mantel, gelijkstroommotor, touchscreenconsole, met laadcellen	SUM1000.9002

Tabel 8.16. Specificaties voor S.U.M.'s van 2.000 I

		Alleen wisselstroommotore	
		Mantel	Zonder mantel
geometrie	Nominaal werkvolume van vloeistof	2.0	00 I
	Minimaal werkvolume van vloeistof	400 I (me mengschroe	t complete efbedekking)
	Totaal kamervolume (vloeistof en gas)	2.7	00
	Diameter BPC-kamer	135 cm	(53 inch)
	Schouderhoogte BPC-kamer	185 cm	(73 inch)
Jgeri	Vloeistofhoogte en nominaal werkvolume	140 cm	(55 inch)
Mei	Vloeistofgeometrie bij verhouding (hoogte/diameter) werkvolume	1	:1
	Ophopingsvolume	<	11
	Algemene mengergeometrie (verhouding hoogte/diameter)	1,1	2:1
	Tankschotten	Ge	en
ef	Mengschroef (hoeveelheid x aantal schoepen)	1:	x 3
chro	Schaal van mengschroef (diameter mengschroef/diameter tank)	1/5	
sbua	Pitch van mengschroefschoepen (hoek)	4!	5°
Ň	Mengschroefdiameter	25,02 cm	(9,85 inch)
	Bereik mengsnelheid	30–35	50 RPM
	Snelheid van tip	39,4–459,7 cm/s ((77,6–904,8 ft/min)
	Mengstroomrichting linksom	Omlaag	pompen
len	Opgeloste hoek roeras	2	7°
hudd	Verschuiving van middellijn roeras	17,8 cm	ו (7 inch)
Sc	Totale lengte aandrijfas	192,5 cm ((75,78 inch)
	Diameter aandrijfas	1,9 cm (0,75 inch)	
	Buitendiameter plastic omhulsel aandrijfas	2,54 cm	ו (1 inch)
	Vrije ruimte mengschroef vanaf tankbodem	5,08 cm	ו (2 inch)
	Motoraandrijving roermechanisme (type, spanning, fase)	Inductie, 208	VAC, 3-fasig
5	Nominaal motorvermogen	745,7 \ Motor: 208	<i>N</i> (1 pk) VAC, 3,2 A
Moto	Nominaal motorkoppel	18 Nm (1	59 in-lb.)
	Tandwielreductie	5	:1
	Methodes voor motorcommunicatie	Op de touchsci communic	reenconsole via atiepoorten

Tabel 8.17. Specificaties voor S.U.M.'s van 2.000 I (vervolg).

		Alleen wisselstroommotoren		
		Mantel	Zonder mantel	
Intel	Manteloppervlak: vol/half volume	5,3 m² (57,3 ft²) / 3,3 m² (35,5 ft²)	-	
ofma	Mantelvolume	26 I (6,9 gal.)	-	
eisto	Manteldebiet bij 3,4 bar (50 psi)	93 l/min (24,5 gal/min)	-	
VIo	Procesverbinding	Hygiënische drieklem 1 inch	-	
regu-	TCU-model: maximale verwarming/koeling	TF24000: 22.500/24.000 W	-	
atuul ring	Geschatte opwarmtijd voor vloeistof (5–37°C)	2,7 uur	-	
ipera le	Geschatte afkoeltijd voor vloeistof (37–5°C)	3,9 uur	-	
Ten	RTD of thermokoppel, 3,18 mm (1/8 inch) OD	Pt-100	(standaard)	
iiner nd)	Totale breedte	180 cm	(70,7 inch)	
ıningsconta ılanken/maı	Totale lengte	171 cm	i (67,4 inch)	
	Algehele hoogte (zonder Powdertainer-arm)	225 cm (88,5 inch)		
ersteı nder p	Gewicht droge skid (massa)	755,1 kg (1664,8 lb.)	551,1 kg (1214,8 lb.)	
0nd (zo	Gewicht natte skid–nominaal werkvolume (massa)	2755,1 kg (6073,8 lb.)	2551,1 kg (5623,8 lb.)	
	Vereiste plafondhoogte voor laden van standaard aandrijfas	292,1 c	m (115 inch)	
	Nominale voeding	100–120 VAC, 50/60 Hz, enkel, 15 A 220–240 VAC, 50/60 Hz, enkel, 10,4 A		
en	Ruisniveau	< 70 dB op 1,5 m		
Algeme	pH- en DO-sonde —autoclaveerbaar type (Applisens, Broadley James, Mettler-Toledo)	12 mm diameter x 215–235 mm invoerlengte x 13,5 PG (pijp) schroefdraad		
	Opslagtemperatuur	-25	5–65°C	
	Relatieve vochtigheid	S.U.M.: 20–80% zonder condensatie Wisselstroommotor: 100% max. (zonder enig dauwcondensatie)		
S	Bereik van bedrijfstemperatuur	2-40 ± 0,1°C (36-104 ± 0,2 °F)		
n ieter:	Motorsnelheid	30–350 RPM		
:vole aram	Volumebereik	500–2.000 I		
anbe iksp:	Maximale statische BPC-druk	0,03 b	ar (0,5 psi)	
A ebruj	Maximale BPC-druk tijdens gebruik	0,007 bar (0,1 psi)		
õ	Continue bedrijfsduur	21 dagen mengen alleen bij nominaal volume		



Afbeelding 8.17. Afmetingen van S.U.M. van 2.000 I (vooraanzicht).



Afbeelding 8.18. Afmetingen van S.U.M. van 2.000 I (bovenaanzicht).

Tabel 8.18. Bestelgegevens voor S.U.M.'s van 2.000 l.

Beschrijving	Cat.nr.
2.000 I zonder mantel, wisselstroommotor, touchscreenconsole, 240 V, met laadcellen	SUM2000.9003
2.000 I met mantel, wisselstroommotor, touchscreenconsole, 240 V, met laadcellen	SUM2000.9004

8.3 Specificaties van de touchscreenconsole

Zie tabel 8.18 hierna voor specificaties van de touchscreenconsole.

		Specificaties		
	Afmetingen (B x H x D)	25 x 52 x 23,4 cm (9,84 x 20,42 x 9,21 inch)		
	Totaal gewicht	18,5 kg (40,8 lb)		
	НМІ	21,34 cm (8,4 inch) lcd-scherm met capacitief touchscreen		
	Constructiemateriaal	AISI 304 roestvrij staal		
en	Elektrische eisen	120/240 VAC, 50/60 Hz, enkel, 15/10,4 A		
geme	Bereik luchtdruk	4,14–6,55 bar (60–95 psi)		
A	Bereik van bedrijfstemperatuur	18–30°C		
	Ruisniveau	< 70 dB op 1,5 m		
	Relatieve vochtigheid	15–80%		
	IP-classificatie	IP54		
	RoHS 2011/65/EU	Conform CE		
	Pompen	Ondersteunt diverse pompen met signaalcontrole van 4-20 mA		
	Laadcellen	Mettler Toledo MTB		
	Ondersteunde pH-sensoren	Thermo, Mettler Toledo, Broadley James		
ita	Ondersteunde geleidbaarheidssensoren	Jumo, Mettler Toledo		
n en da	Knijpventielen	Bimba ACRO 935 knijpventiel 19 mm (3/4 inch) buitendiameter x 3,17 mm (1/8 inch) wandslang		
nente	Alarmen	Ingesteld in de fabriek en door de gebruiker gedefinieerd		
unpor	Communicatiepoorten	USB, Ethernet, Profibus, Modbus RTU		
erne co	E-Stop	Geïntegreerd veiligheidscircuit voor volledig systeem (externe E-Stop ook beschikbaar)		
Exte	Bestandsindelingen	CSV		
	Gegevensregistratie*	Overdracht van door de gebruiker gedefinieerde gegevens via Ethernet, Profibus of Modbus RTU; 72-uurs gegevensopslag voor export via USB		
	Gegevensexport*	Lokaal via USB-flashstation, extern via pc/netwerk met Ethernet, Profibus of Modbus RTU		

*Ontwerp is niet compatibel met CFR11

Tabel 8.20. Meetopties en specificaties van de touchscreenconsole.

	Laadcel	Temperatuur	pН	Geleidbaarheid	BPC en vloeistofdruk in leidingen	Nauwkeurigheid van RPM
Nauwkeurigheid na ± 0,5% van volledige kalibratie schaal		0,2°C	± 0,05 pH	± 5%	± 3,5% van volledige schaal (30 psi)	1 RPM of 1% van instelpunt (grootste van de twee)
Kalibratie	1 tot 3 punten en functie voor nulstellen/tarreren	1 tot 3 punten	1 tot 3 punten	1 tot 3 punten	1 punt; nulstellen/ tarreren	Offset van 1 punt
Resolutie	0,1 kg	0,01°C	0,01 pH	1 µS/cm	0,01 psi.	0,1 RPM
Sensorbereik	50/100 I: 0–300 kg 200 I: 0–600 kg 500 I: 0–900 kg 1.000 I: 0–1500 kg 2.000 I: 0–3300 kg	0–200°C	0–14	20–20.000 µS/cm	0–30 psi	20–350 RPM
Maat- eenheden	kg	Graden Celsius	рН	µS/cm	psi	RPM
Type sonde	3x Mettler Toledo MTB- laadcellen	RTD	Elektrochemisch met S8- connector van 225 mm	Tweepolige geleidbaarheidssensor	Sensor voor eenmalig gebruik, onderdeel van BPC en/of ontwerp met eenheid voor vloeistofoverdracht	N.v.t.

8.4 BPC-specificaties

8.4.1 Standaard BPC's 50 I

Zie de volgende paragrafen voor verschillende tekeningen en specificaties van de standaard S.U.M. BPC van 50 I. De bestelgegevens voor alle typen vindt u in tabel 8.23.

Standaard poeder-op-vloeistof BPC van 50 l (zonder sondepoorten)



Tabel 8.21. BPC van 50 I zonder sondepoorten voor poeder-op-vloeistof toepassingen.

Artikel	Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
1	Afvoer onderkant	12,7 mm (1/2 inch) x 19 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-behuizing met dop van 12,7 mm (1/2 inch)
2	Toevoerleiding	12,7 mm (1/2 inch) x 19 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	Ingestoken 12,7 mm (1/2 inch) MPX-inzetstuk
3	Recirculatie-/ monsternamelijn	12,7 mm (1/2 inch) x 19 mm (3/4 inch) C-Flex slang x 137 cm (54 inch) splitst in 12,7 mm (1/2 inch) x 19 mm (3/4 inch) C-Flex slang x 61 cm (24 inch) en 12,7 mm (1/2 inch) x 19 mm (3/4 inch) C-Flex slang x 30 cm (12 inch)	MPX-invoegstuk met dop van 12,7 mm (1/2 inch), MPX-behuizing met plug van 12,7 mm (1/2 inch)
4	Poedertoevoerpoort	Sanitaire aansluiting 76 mm (3 inch), aansluiting met drie klemmen	Dop met pakking



Standaard vloeistof-op-vloeistof BPC van 50 l (zonder sondepoorten)

Tabel 8.22.	BPC van 50	zonder sondepoorten	vloeistof-op-vloeistof	toepassingen.
-------------	------------	---------------------	------------------------	---------------

Artikel	Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
1	Afvoer onderkant	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-behuizing met dop van 12,7 mm (1/2 inch)
2	Toevoerleiding	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	Ingestoken 12,7 mm (1/2 inch) MPX-inzetstuk
3	Recirculatie-/ monsternamelijn	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 137 cm (54 inch) splitst in 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 61 cm (24 inch) en 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 30 cm (12 inch)	MPX-invoegstuk met dop van 12,7 mm (1/2 inch), MPX- behuizing met plug van 12,7 mm (1/2 inch)
4	Ventilatiefilter	6,4 mm (1/4 inch) x 12,7 mm (1/2 inch) C-Flex-slang x 61 cm (24 inch)	Steriel hydrofoob ventilatiefilter (0,2 micron PVDF—Pall Acro 50)
5	Vulleiding	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	SterilEnz-aansluiting met drie klemmen 38,1 mm (1,5 inch)



Tabel 8.23. BPC van 50 I met sondepoorten voor poeder-op-vloeistof toepassingen.

Artikel	Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
1	Afvoer onderkant	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-behuizing met dop van 12,7 mm (1/2 inch)
2	Toevoerleiding	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	Ingestoken 12,7 mm (1/2 inch) MPX-inzetstuk
3	Recirculatielijn	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 122 cm (48 inch) splitst in 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 61 cm (24 inch) en 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 30 cm (12 inch)	MPX-invoegstuk met dop van 12,7 mm (1/2 inch), MPX-behuizing met plug van 12,7 mm (1/2 inch)
4–5	Sondepoorten (2)	Geen	Aseptische Pall Kleenpak-connectoren KPCHT-serie (vrouwelijk)
6	Thermowell/ monsternamelijn voor kleine volumes	Thermowell-adapter voor diameter van 3,2 mm (1/8 inch) 3,2 mm (1/8 inch) x 6,4 mm (1/4 inch) C-Flex-slang x 30 cm (12 inch)	Luer en SmartSite Kleppoort
7	Poedertoevoerpoort	Sanitaire aansluiting 76 mm (3 inch), aansluiting met drie klemmen	Dop met pakking



Tabal 0.04 DDC van	EO I mot conda	noorton vlooiotof	n vlasistaftas	naaalnaan
Tabel 0.24. DPC van	i bu i met sonde	epoorten vioeistoi-c	p-vioeistoi toe	passingen.

Artikel	Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
1	Afvoer onderkant	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-behuizing met dop van 12,7 mm (1/2 inch)
2	Toevoerleiding	9,5 mm (3/8 inch) x 15,9 mm (5/8 inch) C-Flex-slang x 61 cm (24 inch)	Ingestoken 9,5 mm (3/8 inch) MPX-inzetstuk
3	Recirculatielijn	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch) splitst in 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 61 cm (24 inch) en 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 30 cm (12 inch)	MPX-invoegstuk met dop van 12,7 mm (1/2 inch), MPX-behuizing met plug van 12,7 mm (1/2 inch)
4–5	Sondepoorten (2)	Geen	Aseptische Pall Kleenpak- connectoren KPCHT-serie (vrouwelijk)
6	Thermowell/ monsternamelijn voor kleine volumes	Thermowell-adapter voor diameter van 3,2 mm (1/8 inch) 3,2 mm (1/8 inch) x 6,4 mm (1/4 inch) C-Flex-slang x 30 cm (12 inch)	Luer en SmartSite Kleppoort
7	Ventilatiefilter	6,4 mm (1/4 inch) x 12,7 mm (1/2 inch) C-Flex-slang x 10,2 cm (4 inch)	Dop met pakking
8	Vulleiding	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	SterilEnz-aansluiting met drie klemmen 38,1 mm (1,5 inch)

Standaard vloeistof-op-vloeistof BPC van 50 l (met sondepoorten)

Bestelgegevens voor BPC's van 50 l

Zie tabel 8.23 voor bestelgegevens voor BPC's van 50 l Neem voor meer informatie contact op met uw verkoopvertegenwoordiger.

Opmerking sondepoorten: alle BPC's van 50 l met sondepoorten zijn zo ontworpen dat sondes goed werken bij een turndown van 5:1. Deze BPC's zijn alleen compatibel met de HyPerforma-hardware die in deze publicatie wordt vermeld. Als u een oudere versie van de S.U.M.hardware gebruikt, mag u deze items niet gebruiken. Raadpleeg in plaats daarvan de gebruikershandleiding of gegevensbladen van de verouderde S.U.M.

Tabel 8.25. Bestelgegevens voor BPC's van 50 l.

Afmetingen	Beschrijving	Filmtype	Cat.nr.
50 I	Standaard poeder-op-vloeistof BPC zonder sondepoorten	CX5-14	SH30768.01
50 I	Standaard poeder-op-vloeistof BPC zonder sondepoorten	Aegis5-14	SH30973.01
50	Standaard vloeistof-op-vloeistof BPC zonder sondepoorten	CX5-14	SH30767.01
50	Standaard vloeistof-op-vloeistof BPC zonder sondepoorten	Aegis5-14	SH30983.01
50	Standaard poeder-op-vloeistof BPC met sondepoorten	CX5-14	SH31055.02
50	Standaard poeder-op-vloeistof BPC met sondepoorten	Aegis5-14	SH31051.02
50	Standaard vloeistof-op-vloeistof BPC met sondepoorten	CX5-14	SH31055.04
50	Standaard vloeistof-op-vloeistof BPC met sondepoorten	Aegis5-14	SH31051.01

8.4.2 Standaard BPC's 100 I

Zie de volgende paragrafen voor verschillende tekeningen en specificaties van de standaard S.U.M. BPC van 100 l. De bestelgegevens voor alle typen vindt u in tabel 8.28.

Standaard poeder-op-vloeistof BPC van 100 l (zonder sondepoorten)



Tabel 8.26. BPC van 100 I zonder sondepoorten voor poeder-op-vloeistof toepassingen.

Artikel	Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
1	Afvoer onderkant	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-behuizing met dop van 12,7 mm (1/2 inch)
2	Toevoerleiding	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	Ingestoken 12,7 mm (1/2 inch) MPX-inzetstuk
3	Recirculatie-/ monsternamelijn	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 137 cm (54 inch) splitst in 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 61 cm (24 inch) en 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 30 cm (12 inch)	MPX-invoegstuk met dop van 12,7 mm (1/2 inch), MPX- behuizing met plug van 12,7 mm (1/2 inch)
4	Poedertoevoerpoort	Sanitaire aansluiting 76 mm (3 inch), aansluiting met drie klemmen	Dop met pakking

Standaard vloeistof-op-vloeistof BPC van 100 l (zonder sondepoorten)



Tabel 8.27	. BPC van 100	I zonder sondepoorten	vloeistof-op-vloeistof	toepassingen.
------------	---------------	-----------------------	------------------------	---------------

Artikel	Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
1	Afvoer onderkant	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-behuizing met dop van 12,7 mm (1/2 inch)
2	Toevoerleiding	9,5 mm (3/8 inch) x 15,9 mm (5/8 inch) C-Flex-slang x 61 cm (24 inch)	Ingestoken 9,5 mm (3/8 inch) MPX-inzetstuk
3	Recirculatie-/ monsternamelijn	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 137 cm (54 inch) splitst in 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 61 cm (24 inch) en 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 30 cm (12 inch)	MPX-invoegstuk met dop van 12,7 mm (1/2 inch), MPX- behuizing met plug van 12,7 mm (1/2 inch)
4	Ventilatiefilter	6,4 mm (1/4 inch) x 12,7 mm (1/2 inch) C-Flex-slang x 10,2 cm (4 inch)	Steriel hydrofoob ventilatiefilter (0,2 micron PVDF—Pall Acro 50)
5	Vulleiding	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	SterilEnz-aansluiting met drie klemmen 38,1 mm (1,5 inch)

Standaard poeder-op-vloeistof BPC van 100 l (met sondepoorten)



Tabel 8.28. BPC van 100 l met sondepoorten voor poeder-op-vloeistof toepassingen.

Artikel	Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
1	Afvoer onderkant	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-behuizing met dop van 12,7 mm (1/2 inch)
2	Toevoerleiding	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	Ingestoken 9,5 mm (3/8 inch) MPX-inzetstuk
3	Recirculatielijn	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 137 cm (54 inch) splitst in 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 61 cm (24 inch) en 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 30 cm (12 inch)	MPX-invoegstuk met dop van 12,7 mm (1/2 inch), MPX- behuizing met plug van 12,7 mm (1/2 inch)
4	Thermowell/ monsternamelijn voor kleine volumes	Thermowell-adapter voor diameter van 3,2 mm (1/8 inch) 3,2 mm (1/8 inch) x 6,4 mm (1/4 inch) C-Flex-slang x 30 cm (12 inch)	Luer en SmartSite Kleppoort
5	Poedertoevoerpoort	Sanitaire aansluiting 76 mm (3 inch), aansluiting met drie klemmen	Dop met pakking
6–7	Sondepoorten (2)	Geen	Aseptische Pall Kleenpak- connectoren KPCHT-serie (vrouwelijk)



Standaard vloeistof-op-vloeistof BPC van 100 l

	! f
Tabel 8 29 RPC van 100 i met sondeboorten videistot-ob-vide	istot toenassinden

Artikel	Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
1	Afvoer onderkant	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-behuizing met dop van 12,7 mm (1/2 inch)
2	Toevoerleiding	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	Ingestoken 9,5 mm (3/8 inch) MPX-inzetstuk
3	Recirculatielijn	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 137 cm (54 inch) splitst in 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 61 cm (24 inch) en 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 30 cm (12 inch)	MPX-invoegstuk met dop van 12,7 mm (1/2 inch), MPX- behuizing met plug van 12,7 mm (1/2 inch)
4	Thermowell/ monsternamelijn voor kleine volumes	Thermowell-adapter voor diameter van 3,2 mm (1/8 inch) 3,2 mm (1/8 inch) x 6,4 mm (1/4 inch) C-Flex-slang x 30 cm (12 inch)	Luer en SmartSite Kleppoort
5	Ventilatiefilter	6,4 mm (1/4 inch) x 12,7 mm (1/2 inch) C-Flex-slang x 10,2 cm (4 inch)	Steriel hydrofoob ventilatiefilter (0,2 micron PVDF—Pall Acro 50)
6–7	Sondepoorten (2)	Geen	Aseptische Pall Kleenpak- connectoren KPCHT-serie (vrouwelijk)
8	Vulleiding	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	SterilEnz-aansluiting met drie klemmen 38,1 mm (1,5 inch)

Bestelgegevens voor BPC's van 100 l

Zie tabel 8.28 voor bestelgegevens voor BPC's van 100 l Neem voor meer informatie contact op met uw verkoopvertegenwoordiger.

Opmerking sondepoorten: alle BPC's van 100 I met sondepoorten zijn zo ontworpen dat sondes goed werken bij een turndown van 5:1. Deze BPC's zijn alleen compatibel met de HyPerforma-hardware die in deze publicatie wordt vermeld. Als u een oudere versie van de S.U.M.hardware gebruikt, mag u deze items niet gebruiken. Raadpleeg in plaats daarvan de gebruikershandleiding of gegevensbladen van de verouderde S.U.M.

Tabel 8.30. Bestelgegevens voor BPC's van 100 l.

Afmetingen	Beschrijving	Filmtype	Cat.nr.
100	Standaard poeder-op-vloeistof BPC zonder sondepoorten	CX5-14	SH31046.01
100	Standaard poeder-op-vloeistof BPC zonder sondepoorten	Aegis5-14	SH31052.01
100	Standaard vloeistof-op-vloeistof BPC zonder sondepoorten	CX5-14	SH31046.03
100	Standaard vloeistof-op-vloeistof BPC zonder sondepoorten	Aegis5-14	SH31052.03
100	Standaard poeder-op-vloeistof BPC met sondepoorten	CX5-14	SH31046.02
100	Standaard poeder-op-vloeistof BPC met sondepoorten	Aegis5-14	SH31052.02
100	Standaard vloeistof-op-vloeistof BPC met sondepoorten	CX5-14	SH31046.04
100	Standaard vloeistof-op-vloeistof BPC met sondepoorten	Aegis5-14	SH31052.04

8.4.3 Standaard BPC's 200 I

Zie de volgende paragrafen voor verschillende tekeningen en specificaties van de standaard S.U.M. BPC van 200 I. De bestelgegevens voor alle typen vindt u in tabel 8.33.

Standaard poeder-op-vloeistof BPC van 200 I (zonder sondepoorten)

Voorkant



Tabel 8.31. BPC van 200 l zo	nder sondepoorten voor	r poeder-op-vloeistof	toepassingen.
------------------------------	------------------------	-----------------------	---------------

Artikel	Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
1	Afvoer onderkant	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-inzetstuk met dop van 12,7 mm (1/2 inch)
2	Toevoerleiding	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-behuizing met plug 12,7 mm (1/2 inch)
3	Recirculatie-/ monsternamelijn	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 137 cm (54 inch) splitst in 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 61 cm (24 inch) en 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 30 cm (12 inch)	MPX-invoegstuk met dop van 12,7 mm (1/2 inch), MPX- behuizing met plug van 12,7 mm (1/2 inch)
4	Poedertoevoerpoort	Sanitaire aansluiting 7,6 cm (3 inch), aansluiting met drie klemmen	Dop met pakking



Achterkant

Artikel	Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
1	Afvoer onderkant	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-inzetstuk met dop van 12,7 mm (1/2 inch)
2	Toevoerleiding	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-behuizing met plug 12,7 mm (1/2 inch)
3	Recirculatie-/ monsternamelijn	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 137 cm (54 inch) splitst in 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 61 cm (24 inch) en 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 30 cm (12 inch)	MPX-invoegstuk met dop van 12,7 mm (1/2 inch), MPX-behuizing met plug van 12,7 mm (1/2 inch)
4	Vulleiding	19,1 mm (3/4 inch) x 25,4 mm (1 inch) C-Flex slang x 122 cm (48 inch)	SterilEnz-aansluiting met drie klemmen 38,1 mm (1,5 inch)
5	Ventilatiefilter	6,4 mm (1/4 inch) x 12,7 mm (1/2 inch) C-Flex-slang x 10,2 cm (4 inch)	Steriel hydrofoob ventilatiefilter (0,2 micron PVDF—Pall Acro 50)



Tabel 8.33. BPC van 200 I met sondepoorten voor poeder-op-vloeistof toepassing
--

Artikel	Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
1	Afvoer onderkant	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-behuizing met plug 12,7 mm (1/2 inch)
2	Toevoerleiding	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-behuizing met plug 12,7 mm (1/2 inch)
3	Recirculatie-/ monsternamelijn	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 137 cm (54 inch) splitst in 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 61 cm (24 inch) en 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 30 cm (12 inch)	MPX-invoegstuk met dop van 12,7 mm (1/2 inch), MPX- behuizing met plug van 12,7 mm (1/2 inch)
4	Poedertoevoerpoort	Sanitaire aansluiting 7,6 cm (3 inch), aansluiting met drie klemmen	Dop met pakking
5–6	Sondepoorten (2)	Geen	Aseptische Pall Kleenpak- connectoren KPCHT-serie (vrouwelijk)
7	Thermowell/ monsternamepoort voor kleine volumes	Thermowell-adapter voor diameter van 3,2 mm (1/8 inch) 3,2 mm (1/8 inch) x 6,4 mm (1/4 inch) C-Flex-slang x 30 cm (12 inch)	Luer en SmartSite Kleppoort



Standaard vloeistof-op-vloeistof BPC van 200 l (met sondepoorten)

Tabel 8.34. BPC van	200 I met so	ondepoorten	vloeistof-op-	vloeistof to	epassingen.
	20011110130	macpoonten	viocistor op	1001010110	epussingen.

Artikel	Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
1	Afvoer onderkant	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-inzetstuk met dop van 12,7 mm (1/2 inch)
2	Toevoerleiding	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 6 cm (24 inch)	MPX-behuizing met plug 12,7 mm (1/2 inch)
3	Recirculatie-/ monsternamelijn	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 122 cm (48 inch) splitst in 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 61 cm (24 inch) en 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 30 cm (12 inch)	MPX-invoegstuk met dop van 12,7 mm (1/2 inch), MPX-behuizing met plug van 12,7 mm (1/2 inch)
4	Vulleiding	19,1 mm (3/4 inch) x 25,4 mm (1 inch) C-Flex slang x 122 cm (48 inch)	SterilEnz-aansluiting met drie klemmen 38,1 mm (1,5 inch)
5–6	Sondepoorten (2)	Geen	Aseptische Pall Kleenpak- connectoren KPCHT-serie (vrouwelijk)
7	Thermowell/ monsternamepoort voor kleine volumes	Thermowell-adapter voor diameter van 3,2 mm (1/8 inch) 3,2 mm (1/8 inch) x 6,4 mm (1/4 inch) C-Flex-slang x 30 cm (12 inch)	Luer en SmartSite Kleppoort
8	Ventilatiefilter	6,4 mm (1/4 inch) x 12,7 mm (1/2 inch) C-Flex-slang x 10,2 cm (4 inch)	Steriel hydrofoob ventilatiefilter (0,2 micron PVDF—Pall Acro 50)

Bestelgegevens voor BPC's van 200 l

Zie tabel 8.33 voor bestelgegevens voor BPC's van 200 l Neem voor meer informatie contact op met uw verkoopvertegenwoordiger.

Opmerking sondepoorten: alle BPC's van 200 I met sondepoorten zijn zo ontworpen dat sondes goed werken bij een turndown van 5:1. Deze BPC's zijn alleen compatibel met de HyPerforma-hardware die in deze publicatie wordt vermeld. Als u een oudere versie van de S.U.M.hardware gebruikt, mag u deze items niet gebruiken. Raadpleeg in plaats daarvan de gebruikershandleiding of gegevensbladen van de verouderde S.U.M.

Tabel 8.35. Bestelgegevens voor BPC's van 200 l.

Afmetingen	Beschrijving	Filmtype	Cat.nr.
200	Standaard poeder-op-vloeistof BPC zonder sondepoorten	CX5-14	SH30750.01
200	Standaard poeder-op-vloeistof BPC zonder sondepoorten	Aegis5-14	SH30973.02
200	Standaard vloeistof-op-vloeistof BPC zonder sondepoorten	CX5-14	SH30753.01
200	Standaard vloeistof-op-vloeistof BPC zonder sondepoorten	Aegis5-14	SH30983.02
200	Standaard poeder-op-vloeistof BPC met sondepoorten	CX5-14	SH31054.02
200	Standaard poeder-op-vloeistof BPC met sondepoorten	Aegis5-14	SH31053.02
200	Standaard vloeistof-op-vloeistof BPC met sondepoorten	CX5-14	SH31054.04
200	Standaard vloeistof-op-vloeistof BPC met sondepoorten	Aegis5-14	SH31053.04

8.4.4 Standaard BPC's 500 I

Zie de volgende paragrafen voor verschillende tekeningen en specificaties van de standaard S.U.M. BPC van 500 l. De bestelgegevens voor alle typen vindt u in tabel 8.38.

Standaard poeder-op-vloeistof BPC van 500 l (zonder sondepoorten)



Tabel 8.36. BPC van 500 I zonder sondepoorten voor poeder-op-vloeistof toepassingen.

Artikel	Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
1	Afvoer onderkant	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-inzetstuk met dop van 12,7 mm (1/2 inch)
2	Toevoerleiding	12,7 mm (1/2 inch) x 15,9 mm (5/8 inch) C-Flex-slang x 152 cm (60 inch)	MPX-behuizing met plug 12,7 mm (1/2 inch)
3	Recirculatie-/ monsternamelijn	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 137 cm (54 inch) splitst in 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 61 cm (24 inch) en 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 30 cm (12 inch)	MPX-invoegstuk met dop van 12,7 mm (1/2 inch), MPX-behuizing met plug van 12,7 mm (1/2 inch)
4	Poedertoevoerpoort	Sanitaire aansluiting 7,6 cm (3 inch), aansluiting met drie klemmen	Dop met pakking



Tabel 8.37	. BPC van 500	I zonder sondepoorten	vloeistof-op-vloeistof	toepassingen.
------------	---------------	-----------------------	------------------------	---------------

Artikel	Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
1	Afvoer onderkant	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-inzetstuk met dop van 12,7 mm (1/2 inch)
2	Toevoerleiding	9,5 mm (3/8 inch) x 15,9 mm (5/8 inch) C-Flex-slang x 61 cm (24 inch)	MPX-behuizing met plug 9,5 mm (3/8 inch)
3	Recirculatie-/ monsternamelijn	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 168 cm (66 inch) splitst in 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 61 cm (24 inch) en 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 30 cm (12 inch)	MPX-invoegstuk met dop van 12,7 mm (1/2 inch), MPX-behuizing met plug van 12,7 mm (1/2 inch)
4	Ventilatiefilter	6,4 mm (1/4 inch) x 12,7 mm (1/2 inch) C-Flex-slang x 10,2 cm (4 inch)	Steriel hydrofoob ventilatiefilter (0,2 micron PVDF—Pall Acro 50)
5	Vulleiding	19,1 mm (3/4 inch) x 25,4 mm (1 inch) C-Flex slang x 152 cm (60 inch)	SterilEnz-aansluiting met drie klemmen 38,1 mm (1,5 inch)



Tabel 8.38. BPC van 500 I met sondepoorten voor poeder-op-vloeistof toepassingen.

Artikel	Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
1	Afvoer onderkant	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-behuizing met plug 12,7 mm (1/2 inch)
2	Toevoerleiding	12,7 mm (1/2 inch) x 15,9 mm (5/8 inch) C-Flex-slang x 152 cm (60 inch)	MPX-behuizing met plug 12,7 mm (1/2 inch)
3	Recirculatie-/ monsternamelijn	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 168 cm (66 inch) splitst in 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 61 cm (24 inch) en 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 30 cm (12 inch)	MPX-invoegstuk met dop van 12,7 mm (1/2 inch), MPX-behuizing met plug van 12,7 mm (1/2 inch)
4	Poedertoevoerpoort	Sanitaire aansluiting 7,6 cm (3 inch), aansluiting met drie klemmen	Dop met pakking
5–6	Sondepoorten (2)	Geen	Aseptische Pall Kleenpak- connectoren KPCHT-serie (vrouwelijk)
7	Ongebruikt	Geen	Plug
8	Thermowell/ monsternamepoort voor kleine volumes	Thermowell-adapter voor diameter van 3,2 mm (1/8 inch) 3,2 mm (1/8 inch) x 6,4 mm (1/4 inch) C-Flex-slang x 30 cm (12 inch)	Luer en SmartSite Kleppoort

Standaard poeder-op-vloeistof BPC van 500 I (met sondepoorten)





Tabel 8.39. BPC van 500 I met sondepoorten vloeistof-op-vloeistof toepassingen.

Artikel	Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
1	Afvoer onderkant	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-inzetstuk met dop van 12,7 mm (1/2 inch)
2	Toevoerleiding	9,5 mm (3/8 inch) x 15,9 mm (5/8 inch) C-Flex-slang x 61 cm (24 inch)	MPX-behuizing met plug 9,5 mm (3/8 inch)
3	Recirculatie-/ monsternamelijn	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 168 cm (66 inch) splitst in 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 61 cm (24 inch) en 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 30 cm (12 inch)	MPX-invoegstuk met dop van 12,7 mm (1/2 inch), MPX-behuizing met plug van 12,7 mm (1/2 inch)
4	Ventilatiefilter	6,4 mm (1/4 inch) x 12,7 mm (1/2 inch) C-Flex-slang x 10,2 cm (4 inch)	Steriel hydrofoob ventilatiefilter (0,2 micron PVDF—Pall Acro 50)
5	Vulleiding	19,1 mm (3/4 inch) x 25,4 mm (1 inch) C-Flex slang x 152 cm (60 inch)	SterilEnz-aansluiting met drie klemmen 38,1 mm (1,5 inch)
6–7	Sondepoorten (2)	Geen	Aseptische Pall Kleenpak- connectoren KPCHT-serie (vrouwelijk)
8	Dompelhuls/ monster met klein volume	Thermowell-adapter voor diameter van 3,2 mm (1/8 inch) 3,2 mm (1/8 inch) x 6,4 mm (1/4 inch) C-Flex-slang x 30 cm (12 inch)	Luer en SmartSite Kleppoort

Thermo Scientific

Bestelgegevens voor BPC's van 500 l

Zie onderstaande tabel 8.40 voor bestelgegevens voor BPC's van 500 I Neem voor meer informatie contact op met uw verkoopvertegenwoordiger.

Tabel 8.40. Bestelgegevens voor BPC's van 500 l.

Afmetingen	Beschrijving	Filmtype	Cat.nr.
500 I	Standaard poeder-op-vloeistof BPC zonder sondepoorten	CX5-14	SH30751.01
500 I	Standaard poeder-op-vloeistof BPC zonder sondepoorten	Aegis5-14	SH30973.03
500 I	Standaard vloeistof-op-vloeistof BPC zonder sondepoorten	CX5-14	SH30754.01
500 I	Standaard vloeistof-op-vloeistof BPC zonder sondepoorten	Aegis5-14	SH30983.03
500	Standaard poeder-op-vloeistof BPC met sondepoorten	CX5-14	SH30751.02
500 I	Standaard poeder-op-vloeistof BPC met sondepoorten	Aegis5-14	SH30974.03
500	Standaard vloeistof-op-vloeistof BPC met sondepoorten	CX5-14	SH30754.02
500	Standaard vloeistof-op-vloeistof BPC met sondepoorten	Aegis5-14	SH30982.03

8.4.5 Standaard BPC's 1.000 I

Zie de volgende paragrafen voor verschillende tekeningen en specificaties van de standaard S.U.M. BPC van 1.000 l. De bestelgegevens voor alle typen vindt u in tabel 8.43.





Tabel 8.41. BPC van 1.000 I zonder sondepoorten voor poeder-op-vloeistof toepassingen.

Artikel	Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
1	Afvoer onderkant	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-inzetstuk met dop van 12,7 mm (1/2 inch)
2	Toevoerleiding	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 183 cm (72 inch)	MPX-behuizing met plug 12,7 mm (1/2 inch)
3	Recirculatie-/ monsternamelijn	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 198 cm (78 inch) splitst in 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 61 cm (24 inch) en 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 30,5 cm (12 inch)	MPX-invoegstuk met dop van 12,7 mm (1/2 inch), MPX-behuizing met plug van 12,7 mm (1/2 inch)
4	Poedertoevoerpoort	Sanitaire aansluiting 7,6 cm (3 inch), aansluiting met drie klemmen	Dop met pakking

Standaard vloeistof-op-vloeistof BPC van 1.000 l (zonder sondepoorten)



Achterkant

Tabel 8.42. BPC van 1.000 I zonder sondepoorten vloeistof-op-vloeistof toepassingen.

Artikel	Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
1	Afvoer onderkant	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-inzetstuk met dop van 12,7 mm (1/2 inch)
2	Toevoerleiding	9,5 mm (3/8 inch) x 15,9 mm (5/8 inch) C-Flex-slang x 61 cm (24 inch)	MPX-behuizing met plug 9,5 mm (3/8 inch)
3	Recirculatie-/ monsternamelijn	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 198 cm (78 inch) splitst in 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 61 cm (24 inch) en 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 30 cm (12 inch)	MPX-invoegstuk met dop van 12,7 mm (1/2 inch), MPX-behuizing met plug van 12,7 mm (1/2 inch)
4	Ventilatiefilter	6,4 mm (1/4 inch) x 12,7 mm (1/2 inch) C-Flex-slang x 10,2 cm (4 inch)	Steriel hydrofoob ventilatiefilter (0,2 micron PVDF—Pall Acro 50)
5	Vulleiding	19,1 mm (3/4 inch) x 25,4 mm (1 inch) C-Flex slang x 183 cm (72 inch)	SterilEnz-aansluiting met drie klemmen 38,1 mm (1,5 inch)

Standaard poeder-op-vloeistof BPC van 1.000 l (met sondepoorten)



Tabel 8.43. BPC van 1.000 I met sondepoorten voor poeder-op-vloeistof toepassingen.

Artikel	Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
1	Afvoer onderkant	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-behuizing met plug 12,7 mm (1/2 inch)
2	Toevoerleiding	12,7 mm (1/2 inch) x 15,9 mm (5/8 inch) C-Flex-slang x 122 cm (48 inch)	MPX-behuizing met plug 12,7 mm (1/2 inch)
3	Recirculatie-/ monsternamelijn	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 198 cm (78 inch) splitst in 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 61 cm (24 inch) en 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 30 cm (12 inch)	MPX-invoegstuk met dop van 12,7 mm (1/2 inch), MPX-behuizing met plug van 12,7 mm (1/2 inch)
4	Poedertoevoerpoort	Sanitaire aansluiting 7,6 cm (3 inch), aansluiting met drie klemmen	Dop met pakking
5–6	Sondepoorten (2)	Geen	Aseptische Pall Kleenpak- connectoren KPCHT-serie (vrouwelijk)
7	Ongebruikt	Geen	Plug
8	Thermowell/ monsternamepoort voor kleine volumes	Thermowell-adapter voor diameter van 3,2 mm (1/8 inch) 3,2 mm (1/8 inch) x 6,4 mm (1/4 inch) C-Flex-slang x 30 cm (12 inch)	Luer en SmartSite Kleppoort

Standaard vloeistof-op-vloeistof BPC van 1.000 l (met sondepoorten)





Tabel 8.44. BPC van 1.000 I met sondepoorten vloeistof-op-vloeistof toepassingen.

Artikel	Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
1	Afvoer onderkant	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-inzetstuk met dop van 12,7 mm (1/2 inch)
2	Toevoerleiding	9,5 mm (3/8 inch) x 15,9 mm (5/8 inch) C-Flex-slang x 61 cm (24 inch)	MPX-behuizing met plug 9,5 mm (3/8 inch)
3	Recirculatie-/ monsternamelijn	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 198 cm (78 inch) splitst in 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 61 cm (24 inch) en 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 30 cm (12 inch)	MPX-invoegstuk met dop van 12,7 mm (1/2 inch), MPX-behuizing met plug van 12,7 mm (1/2 inch)
4	Ventilatiefilter	6,4 mm (1/4 inch) x 12,7 mm (1/2 inch) C-Flex-slang x 10,2 cm (4 inch)	Steriel hydrofoob ventilatiefilter (0,2 micron PVDF—Pall Acro 50)
5	Vulleiding	19,1 mm (3/4 inch) x 25,4 mm (1 inch) C-Flex slang x 183 cm (72 inch)	SterilEnz-aansluiting met drie klemmen 38,1 mm (1,5 inch)
6–7	Sondepoorten (2)	Geen	Aseptische Pall Kleenpak- connectoren KPCHT-serie (vrouwelijk)
8	Ongebruikt	Geen	Plug
9	Dompelhuls/ monster met klein volume	Thermowell-adapter voor diameter van 3,2 mm (1/8 inch) 3,2 mm (1/8 inch) x 6,4 mm (1/4 inch) C-Flex x 30 cm (12 inch)	Luer en SmartSite Kleppoort

Bestelgegevens voor BPC's van 1.000 l

Zie onderstaande tabel 8.43 voor bestelgegevens voor BPC's van 1.000 I Neem voor meer informatie contact op met uw verkoopvertegenwoordiger.

Tabel 8.45. Bestelgegevens voor BPC's van 1.000 l.

Afmetingen	Beschrijving	Filmtype	Cat.nr.
1.000 I	Standaard poeder-op-vloeistof BPC zonder sondepoorten	CX5-14	SH30752.02
1.000 l	Standaard poeder-op-vloeistof BPC zonder sondepoorten	Aegis5-14	SH30973.04
1.000 I	Standaard vloeistof-op-vloeistof BPC zonder sondepoorten	CX5-14	SH30755.01
1.000 I	Standaard vloeistof-op-vloeistof BPC zonder sondepoorten	Aegis5-14	SH30983.04
1.000 I	Standaard poeder-op-vloeistof BPC met sondepoorten	CX5-14	SH30752.02
1.000 I	Standaard poeder-op-vloeistof BPC met sondepoorten	Aegis5-14	SH30974.04
1.000 I	Standaard vloeistof-op-vloeistof BPC met sondepoorten	CX5-14	SH30755.02
1.000 I	Standaard vloeistof-op-vloeistof BPC met sondepoorten	Aegis5-14	SH30982.04

8.4.6 Standaard BPC's 2.000 I

Zie de volgende paragrafen voor verschillende tekeningen en specificaties van de standaard S.U.M. BPC van 2.000 I. De bestelgegevens voor alle typen vindt u in tabel 8.48.



Poeder-op-vloeistof BPC van 2.000 l (zonder sondepoorten)

Tabel 8.46. BPC van 2.000 I zonder sondepoorten voor poeder-op-vloeistof toepassingen.

Artikel	Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
1	Afvoer onderkant	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-inzetstuk met dop van 12,7 mm (1/2 inch)
2	Toevoerleiding	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 183 cm (72 inch)	MPX-behuizing met plug 12,7 mm (1/2 inch)
3	Recirculatie-/ monsternamelijn	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 198 cm (78 inch) splitst in 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 61 cm (24 inch) en 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 30,5 cm (12 inch)	MPX-invoegstuk met dop van 12,7 mm (1/2 inch), MPX-behuizing met plug van 12,7 mm (1/2 inch)
4	Poedertoevoerpoort	Sanitaire aansluiting 7,6 cm (3 inch), aansluiting met drie klemmen	Dop met pakking

Standaard vloeistof-op-vloeistof BPC van 2.000 l (zonder sondepoorten)



Tabel 8.47	. BPC van 2.000	I zonder sondepoorten	vloeistof-op-vloeistof	toepassingen.
------------	-----------------	-----------------------	------------------------	---------------

Artikel	Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
1	Afvoer onderkant	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-inzetstuk met dop van 12,7 mm (1/2 inch)
2	Toevoerleiding	9,5 mm (3/8 inch) x 15,9 mm (5/8 inch) C-Flex-slang x 61 cm (24 inch)	MPX-behuizing met plug 9,5 mm (3/8 inch)
3	Recirculatie-/ monsternamelijn	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 198 cm (78 inch) splitst in 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 61 cm (24 inch) en 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 30 cm (12 inch)	MPX-invoegstuk met dop van 12,7 mm (1/2 inch), MPX-behuizing met plug van 12,7 mm (1/2 inch)
4	Ventilatiefilter	6,4 mm (1/4 inch) x 12,7 mm (1/2 inch) C-Flex-slang x 10,2 cm (4 inch)	Steriel hydrofoob ventilatiefilter (0,2 micron PVDF—Pall Acro 50)
5	Vulleiding	19,1 mm (3/4 inch) x 25,4 mm (1 inch) C-Flex slang x 183 cm (72 inch)	SterilEnz-aansluiting met drie klemmen 38,1 mm (1,5 inch)



Standaard poeder-op-vloeistof BPC van 2.000 l (met sondepoorten)

Tabel 8.48. BPC van 2.000 I met sondepoorten voor poeder-op-vloeistof toepassingen.

Artikel	Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
1	Afvoer onderkant	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-behuizing met plug 12,7 mm (1/2 inch)
2	Toevoerleiding	12,7 mm (1/2 inch) x 15,9 mm (5/8 inch) C-Flex-slang x 122 cm (48 inch)	MPX-behuizing met plug 12,7 mm (1/2 inch)
3	Recirculatie-/ monsternamelijn	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 198 cm (78 inch) splitst in 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 61 cm (24 inch) en 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 30 cm (12 inch)	MPX-invoegstuk met dop van 12,7 mm (1/2 inch), MPX-behuizing met plug van 12,7 mm (1/2 inch)
4	Poedertoevoerpoort	Sanitaire aansluiting 7,6 cm (3 inch), aansluiting met drie klemmen	Dop met pakking
5–6	Sondepoorten (2)	Geen	Aseptische Pall Kleenpak- connectoren KPCHT-serie (vrouwelijk)
7	Ongebruikt	Geen	Plug
8	Thermowell/ monsternamepoort voor kleine volumes	Thermowell-adapter voor diameter van 3,2 mm (1/8 inch) 3,2 mm (1/8 inch) x 6,4 mm (1/4 inch) C-Flex-slang x 30 cm (12 inch)	Luer en SmartSite Kleppoort

Standaard vloeistof-op-vloeistof BPC van 2.000 l (met sondepoorten)

Voorkant



Tabel 8.49. BPC van 2.000 I met sondepoorten vloeistof-op-vloeistof toepassingen.

Artikel	Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
1	Afvoer onderkant	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex-slang x 122 cm (48 inch)	MPX-inzetstuk met dop van 12,7 mm (1/2 inch)
2	Toevoerleiding	9,5 mm (3/8 inch) x 15,9 mm (5/8 inch) C-Flex-slang x 61 cm (24 inch)	MPX-behuizing met plug 9,5 mm (3/8 inch)
3	Recirculatie-/ monsternamelijn	12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 198 cm (78 inch) splitst in 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 61 cm (24 inch) en 12,7 mm (1/2 inch) x 19,1 mm (3/4 inch) C-Flex slang x 30 cm (12 inch)	MPX-invoegstuk met dop van 12,7 mm (1/2 inch), MPX-behuizing met plug van 12,7 mm (1/2 inch)
4	Ventilatiefilter	6,4 mm (1/4 inch) x 12,7 mm (1/2 inch) C-Flex-slang x 10,2 cm (4 inch)	Steriel hydrofoob ventilatiefilter (0,2 micron PVDF—Pall Acro 50)
5	Vulleiding	19,1 mm (3/4 inch) x 25,4 mm (1 inch) C-Flex slang x 183 cm (72 inch)	SterilEnz-aansluiting met drie klemmen 38,1 mm (1,5 inch)
6–7	Sondepoorten (2)	Geen	Aseptische Pall Kleenpak- connectoren KPCHT-serie (vrouwelijk)
8	Ongebruikt	Geen	Plug
9	Dompelhuls/ monster met klein volume	Thermowell-adapter voor diameter van 3,2 mm (1/8 inch) 3,2 mm (1/8 inch) x 6,4 mm (1/4 inch) C-Flex-slang x 30 cm (12 inch)	Luer en SmartSite Kleppoort
Bestelgegevens voor BPC's van 2.000 I

Zie onderstaande tabel 8.43 voor bestelgegevens voor BPC's van 2.000 I Neem voor meer informatie contact op met uw verkoopvertegenwoordiger.

Tabel 8.50. Bestelgegevens voor BPC's van 2.000 I.

Afmetingen	Beschrijving	Filmtype	Cat.nr.
2.000 I	Standaard poeder-op-vloeistof BPC zonder sondepoorten	CX5-14	SH30770.01
2.000 I	Standaard poeder-op-vloeistof BPC zonder sondepoorten	Aegis5-14	SH30973.05
2.000	Standaard vloeistof-op-vloeistof BPC zonder sondepoorten	CX5-14	SH30769.01
2.000	Standaard vloeistof-op-vloeistof BPC zonder sondepoorten	Aegis5-14	SH30983.05
2.000 I	Standaard poeder-op-vloeistof BPC met sondepoorten	CX5-14	SH30770.02
2.000 I	Standaard poeder-op-vloeistof BPC met sondepoorten	Aegis5-14	SH30974.05
2.000 I	Standaard vloeistof-op-vloeistof BPC met sondepoorten	CX5-14	SH30769.02
2.000	Standaard vloeistof-op-vloeistof BPC met sondepoorten	Aegis5-14	SH30982.05

8.4.7 Standaard open-top liners en mengschroefhulzen

Zie de onderstaande tekeningen en tabellen 8.49–8.51 voor specificaties en bestelgegevens voor open-top liners en mengschroefhulzen.



Afvoer onderkant

Tabel 8.51. Standaard open-top liner.

Beschrijving	Buizenset (ID x OD x lengte)	Eindbehandeling
Afvoer onderkant	182,9 mm (72 inch) van 12,7 mm (1/2 inch) C-Flex slangen	Connector voor MPX- invoegstuk



Let op: de lagerpot voor open-top mengen wordt standaard samen met de tankhardware geleverd. De eindbehandeling voor de mengschroef voor 2.000 l heeft een aansluiting met drie klemmen in plaats van een snelaansluiting.

Bestelgegevens voor open-top liners en mengschroefhulzen

Afmetingen	Beschrijving	Filmtype	Cat.nr.
50	Standaard open-top liner	CX3-9	SH30762.04
100 I	Standaard open-top liner	CX3-9	SH30762.06
200	Standaard open-top liner	CX3-9	SH30762.01
500 I	Standaard open-top liner	CX3-9	SH30762.02
1.000 I	Standaard open-top liner	CX3-9	SH30762.03
2.000 I	Standaard open-top liner	CX3-9	SH30762.05

Tabel 8.52. Bestelgegevens voor open-top liners voor de S.U.M.

	Tabel 8.53. Bestelgegevens	voor mengschroefhulzer	ו voor de S.U.M.
--	----------------------------	------------------------	------------------

Afmetingen	Beschrijving	Cat.nr.
50	Mengschroefhuls voor open-top mengen	SH30749.06
100	Mengschroefhuls voor open-top mengen	SH30749.14
200	Mengschroefhuls voor open-top mengen	SH30749.08
500 I	Mengschroefhuls voor open-top mengen	SH30749.10
1.000	Mengschroefhuls voor open-top mengen	SH30749.10
2.000	Mengschroefhuls voor open-top mengen	SH30772.01

8.4.8 Aangepaste BPC-producten

Zie tabel 8.52 hierna voor beschrijvingen van verschillende aangepaste BPC-producten. Neem voor meer informatie contact op met uw verkoopvertegenwoordiger.

Let op: Niet alle opties zijn beschikbaar voor alle poorten. Poorttype, poortlocatie, kamerafmetingen en mengeenheid zijn niet aanpasbaar. Zie voor meer informatie de selectiegidsen in de BPC-catalogus.

 Tabel 8.54. Specificatiegegevens voor aangepaste BPC-producten.

Categorie	Opties/capaciteiten	Opmerkingen
Buistype	C-Flex, platinum-cured silicone, PVC, PharMed, PharmaPure	Meer informatie is beschikbaar in de Keuzegids voor buizen.
Grootte van de buis	Bereiken met binnendiameters van 3,18–25,4 mm (1/8–1 inch) in verschillende lengtes	Meer informatie is beschikbaar in de Keuzegids voor buizen.
Connectoren	Luers, CPC-snelaansluitingen, SIP-connectoren, drieklem, Kleenpak, Lynx, SmartSite, Clave, Lynx steam thru, CPC steam thru, Gore stoomklep, Gore Mini TC, BioQuate, SterilEnz, eindplug enz.	Meer informatie is beschikbaar in de Keuzegids voor aansluitsystemen.
Sondepoorten/ lijntoevoerpoorten	Poorten kunnen worden toegevoegd als ze compatibel zijn met de hardware.	De herbruikbare sondeconnector is geschikt voor Kleenpak- of AseptiQuik-connectoren
Wegwerpsensoren	Druksensor: PendoTECH- en Thermo Scientific- oplossingen (PendoTECH is standaard aanwezig op 500 l en 1.000 l) DO en pH: Thermo Scientific en PreSens pH: Mettler Toledo	Keuze aan gekwalificeerde leveranciers is beschikbaar
Poortmaten	Enkel beperkte engineer-to-order aanpassing	Afhankelijk van de locatie in de BPC en pasvorm binnen de hardware (bijv. een poort met een diameter van 2,54 cm (1 inch) op een aftapleiding)
Herschikking van lijnen aan bestaande poorten	Beperkte aanpassingsmogelijkheden, zoals verplaatsing van de monstername-/ thermowellpoort van een sondeslangpoort, of omwisseling van uitlaatleiding en vloeistofleidingen	Afhankelijk van de locatie in de BPC en pasvorm binnen de hardware
Lijnen met ondergedompelde buizen	Beperkte aanpassing mogelijk	Lengte kan de mengschroef en as niet storen
Filters aan media- en supplementinlaten	Alleen beperkte engineer-to-order aanpassing	N.v.t.

Buitenverpakking	'Plat verpakt' geleverd met twee polyethyleen buitenlagen	
Label	Beschrijving, productcode, partijnummer, vervaldatum op buitenverpakking, verzenddoos	
Gammastraling	Straling (25 tot 40 kGy) in buitenverpakking	
Verzonden container	Duurzame kartonnen doos	
Documentatie	Analysecertificaat wordt voor elke levering bij elke partij verschaft	

Tabel 8.55.	Informatie	over BPC	-verpakking

8.5 Specificaties van accessoires en opties

De volgende paragrafen bevatten informatie over accessoires en opties voor de S.U.M.

8.5.1 Laadcellen

Laadcellen werken als een mechanisch ophang- en weegsysteem op de S.U.M. Laadcellen kunnen worden aangeschaft op het moment van verkoop van de apparatuur of als retro-fit kits voor bestaande S.U.M.-eenheden. Laadcelsystemen bestaan uit drie laadcellen, een optelblok, een display en bedrading.

Laadcellen uit de Mettler Toledo 0745A-serie worden gebruikt om met S.U.M.-eenheden van 2.000 I te kunnen wegen (afbeelding 8.19). Mettler Toledo MTB-laadcellen (afbeelding 8.20 en 8.21) worden gebruikt om met alle andere S.U.M.-eenheden te kunnen wegen.



Afbeelding 8.19. Mettler Toledo 0745A-laadcel voor de S.U.M. van 2.000 l





Afbeelding 8.20. Mettler Toledo MTBlaadcel voor S.U.M-eenheden van 50-1.000 I

Afbeelding 8.21. Close-up van Mettler Toledo MTB-laadcel.

Laadcellen worden gewoonlijk radiaal gemonteerd in sets van drie. De montagelocatie is iets anders voor elke grootte om de onderste afvoer- of spoelsystemen en -slangen gemakkelijker toegankelijk te maken. Zie tabel 8.54 voor specificaties van de laadcellen en tabel 8.55 voor bestelgegevens.

Tabel 8.56. Specificatiegegevens voor S.U.M.-laadcellen.

Parameter weegmodule	S.U.M. 50–100 I	S.U.M. 200 I	S.U.M. 500 I	S.U.M. 1.000 I	S.U.M. 2.000 I
Modelnummer	MTB-100	MTB-200	MTB-300	MTB-500	0745A
Nominale capaciteit per laadcel	100 kg (220 lb) 200 kg (441 lb) 300 kg (661 lb) 500 kg (1102 lb)		1100 kg (2500 lb)		
Veilige laadlimiet	150% van nominale capaciteit				
Veilige laadlimiet zijkant	100% van nominale capaciteit				
Veilige dynamische lading	70% van nominale capaciteit				
Gewicht (inclusief laadcel),	0.6 kg (1.3 lb)		0.6 kg (1.3 lb)		Volledige eenheid 7,2 kg (16 lb)
nominaal		Alleen laadcel 0,9 kg (1,98 lb)			
Materiaal	304 roestvrij staal				

Afmetingen	Beschrijving	Cat.nr.
50, 100 l	(3) Laadcel met kabels	SV51145.01
200 I	(3) Laadcel met kabels	SV51145.02
500 I	(3) Laadcel met kabels	SV51145.03
1.000	(3) Laadcel met kabels	SV51145.05
2.000	(3) Laadcel met kabels	SV51146.04

Tabel 8.57. Bestelgegevens voor laadcelkits van Mettler Toledo.

8.5.2 Powdertainer-arm

De Powdertainer-arm is als optie beschikbaar voor poeder-opvloeistof toepassingen. Powdertainer-armen zijn verkrijgbaar in twee maten: één voor eenheden van 50 tot 1.000 l, en één voor eenheden van 2.000 l (afbeelding 8.22). De arm houdt de poedercontainer boven de mixer vast en bevestigt deze met een klem aan de BPC. De arm kan verticaal worden aangepast en is draaibaar zodat de poedercontainer handig op de hanger kan worden getild. Zie tabel 8.56 voor bestelgegevens voor de Powdertainer-arm.



Afbeelding 8.22. Powdertainer-armen voor eenheden van 50–1.000 I (links) en eenheden van 2.000 I (rechts)

Afmetingen	Beschrijving	Cat.nr.
50–1.000 I	Powdertainer-arm voor mixers van 50-1.000 I	SV51002.01
2.000 I	Powdertainer-arm voor mixers van 2.000 l	SV51002.02

Tabel 8.58. Bestelgegevens voor Powdertainer-armen.

8.5.3 Kabelbeheersysteem

Het kabelbeheersysteem (afbeelding 8.23) is een accessoire dat gebruikers helpt de draden en kabels uit de touchscreenconsole bij elkaar te houden. Indien bij uw systeem besteld, zijn de kabelbeheerhaken bij levering al aan de mixer bevestigd. Kabelbeheersystemen kunnen ook achteraf aan bestaande S.U.M.eenheden worden toegevoegd. Zie tabel 8.57 voor bestelgegevens voor kabelbeheersystemen.



Afbeelding 8.23. Kabelbeheerarm.

Tabel 8.59. Bestelgegevens voor kabelbeheersysteem.

Afmetingen	Beschrijving	Cat.nr.
50–2.000 I	Kabelbeheerarm	SV50992.12

8.5.4 Diverse items

De volgende hulpcomponenten ondersteunen de werking van de HyPerforma S.U.M. en verbeteren de prestaties van het hele systeem. Bestelgegevens voor alle verschillende items vindt u in tabel 8.58.

Autoclaaftray voor sondes

De autoclaafplaat (afbeelding 8.24) bevat de elektrochemische sondes en balg die worden geplaatst tijdens het autoclaaf-sterilisatieproces. De tray biedt ruimte voor twee sondes.

Ontwerpelementen zijn:

- Gemaakt van roestvrij staal
- Een plastic handgreep voor gemakkelijk transport meteen vanuit de autoclaaf
- Mogelijkheid om sondes onder een hoek van 15% te plaatsen voor een grotere levensduur van sonde/membraan
- Mogelijkheid om te voorkomen dat de sondebalgen samenvouwen tijdens de sterilisatie



Afbeelding 8.24. Autoclaaftray voor aseptische Kleenpakconnectoren.

Robuuste slangklemmen

Robuuste slangklemmen (afbeelding 8.25) worden gebruikt om ongebruikte lijnsets dicht te knijpen en te voorkomen dat procesvloeistoffen ontsnappen. Slangklemmen moeten op hun plek zitten voordat steriele sondes worden ingebracht, zodat de sondepoorten afgesloten zijn.



Afbeelding 8.25. Slangklem.

Knijpventielen

Knijpventielen (afbeelding 8.26) worden aangesloten op de touchscreenconsole en worden gebruikt om de vloeistofstroom door de BPC-slangen automatisch te reguleren tijdens het vullen of aftappen.



Afbeelding 8.26. Knijpventiel.

Sondeclips

Plastic sondeclips (afbeelding 8.27) worden gebruikt om sondes op de S.U.M.-tank op hun plek te houden. De sondeclips kunnen onafhankelijk worden verplaatst en zijn aangebracht op een dunne beugel boven de uitsparing in de tank voor sondepoorten. Plastic sondeclips worden bevestigd door ze met stevige druk op de beugel aan te brengen.



Afbeelding 8.27. Plastic sondeclip.

Pompen

Pompen zijn optioneel voor S.U.M.-systemen. Indien gebruikt in combinatie met de touchscreenconsole, stellen pompen gebruikers in staat vloeistof automatisch bij te vullen en af te tappen en bolussen tijdens het mengen toe te voeren. Verschillende pompmaten zijn verkrijgbaar om in te spelen op de verschillende behoeften van gebruikers (afbeelding 8.28). Grote pompen worden gebruikt om vloeistof uit de S.U.M. aan te vullen en af te tappen. Kleine pompen worden hoofdzakelijk gebruikt om bufferoplossingen en pH-bolussen tijdens het mengen toe te dienen.



Afbeelding 8.28. Watson-Marlowpompen op planken.

Planken

Planken om pompen op te bergen (afbeelding 8.29) zijn een optioneel accessoire voor S.U.M.-systemen. Er zijn drie schappen gemonteerd op de arm die op de tank is bevestigd. Ze kunnen één grote en twee kleine Watson-Marlow-pompen dragen.



Mand

Een mand is als accessoire beschikbaar voor de S.U.M. De mand biedt opslagruimte voor flessen met buffer- of pH-oplossingen (afbeelding 8.30). De mand wordt gemonteerd op de arm die boven de pompplanken is bevestigd.



Afbeelding 8.30. Mand voor oplossingflessen.

Bestelgegevens

Zie tabel 8.58 voor bestelgegevens voor diverse items.

Tabel 8.60. Bestelgegevens voor diverse items.

Beschrijving	Cat.nr.
Autoclaaftray voor aseptische connectoren.	SV50177.01
Sonde-eenheid (niet-steriel, voor gebruik in autoclaaf)	SH30720.01
Sondeclips (4)	SV50177P.01
Robuuste slangklemmen (enkel)	SV20664.01
Robuuste slangklemmen (10)	SV20664.04
Watson-Marlow 120 V wisselstroompomp — groot (0,004–6,85 l/min.)	SV511054.01
Watson-Marlow 240 V wisselstroompomp — groot (0,004–6,85 l/min.)	SV511054.02
Watson-Marlow 120 – 240 wisselstroompomp — klein (0,09–170 ml/min.)	SV511054.03
Pompplanken en mand	SV50236A.12
T-stekker — RS485/Modbus/Canbus	SV51142.901
Terminator — RS485/Modbus/Canbus	SV51142.900
T-stekker — Profibus	SV51142.902
Terminator — Profibus	SV51142.903
Eindstekker, externe E-Stop	SV51142.904
pH-sonde (Thermo Scientific), 225 mm K8	SV51147.02
pH-sonde (Broadley James), 225 mm K8	SV51147.03
pH-sonde (Mettler Toledo), 225 mm K8	SV51147.01
RTD met dubbel element	SV50999.08
Geleidbaarheidssonde (Jumo)	SV51148.01
Geleidbaarheidssonde (Mettler Toledo)	SV51148.02
Knijpventiel, vulleiding	SV51108.05
Knijpventiel, aftapleiding	SV51108.08
Kabel, hulpingang/uitgang 4-20 mA	SV51142.01
Kabel, externe E-Stop	SV51142.02
Kabel, extern alarm	SV51142.03
Kabel, verouderd 17-pins	SV51142.04
Kabel, Modbus TCP	SV51142.100
Kabel, E-Box naar TCU (Lauda)	SV51142.101
Kabel, E-Box naar TCU (Neslab)	SV51142.102
Kabel, Profibus	SV51142.103
Kabel, RS485/Modbus RTU	SV51142.104
Kabel, CANBus	SV51142.106
Kabel, enkele RTD	SV51142.200
Kabel, dubbele RTD Lemo-connector (voor Lauda TCU)	SV51142.201
Kabel, dubbele RTD DB9-connector (voor Neslab TCU)	SV51142.202
Kabel, druksensor	SV51142.300
Kabel, pH voor K8S-connector	SV51142.400
Kabel, pH voor VP6-connector	SV51142.401
Kabel, geleidbaarheidssensor met VarioPin-connector (Mettler Toledo)	SV51142.500
Kabel, geleidbaarheidssensor met M12-connector (Jumo)	SV51142.501

Beschrijving	Cat.nr.
Kabel, gelijkstroommotor D-Sub, comm, en vermogen	SV51142.600
Kabel, hulpuitgang voor wisselstroom	SV51142.800
Kabel, voedingskabel voor wisselstroommotor	SV50986.18
Motor, gelijkstroom met geïntegreerde aandrijving, 400 W, overbrengingsverhouding 7,5:1	SV51152.01
Motor, wisselstroom, 1 pk, overbrengingsverhouding 5:1	SV50237.29
Touchscreenconsole	SV51151.01

Tabel 8.61. Bestelgegevens voor diverse items (vervolg).

Reserveonderdelen

Zie de bij het ETP geleverde tekeningen voor informatie over aanbevolen reserveonderdelen.

8.6 Configureerbare opties

Zie afbeelding 8.31–8.37 hierna voor configureerbare opties voor de HyPerforma S.U.M. met touchscreenconsole.



Afbeelding 8.31. Verplicht configureerbare opties voor S.U.M.



Afbeelding 8.32. Motoropties.







Afbeelding 8.34. Opties voor pH-sensor.



Afbeelding 8.35. Opties voor geleidbaarheidssensor.



Afbeelding 8.36. Opties voor voedingskabel.





Afbeelding 8.38. Pompopties.



Afbeelding 8.39. Opties voor secundaire montage.



Afbeelding 8.40. Opties voor knijpventielen.

0

Algemene bestelinformatie

Inhoudsopgave

- 9.1 Bestelinstructies
- 9.2 Contactgegevens voor bestellen en ondersteuning
- 9.3 Technische ondersteuning

9.1 Bestelinstructies

BPC's en hardwarecomponenten voor de HyPerforma S.U.M. kunnen rechtstreeks besteld worden bij Thermo Fisher Scientific. Deze artikelen bestaan uit alle componenten met onderdeelnummers die beginnen met de volgende tekens:

- SH
- SV
- SUM

9.2 Contactgegevens voor bestellen en ondersteuning

In Noord-, Midden- Zuid-Amerika en Azië

1726 Hyclone Drive Logan, Utah 84321 Verenigde Staten Tel: +1 435 792 8500 E-mail: customerservice.bioprocessing@thermofisher.com

In Europa

Unit 9 Atley Way Cramlington, NE 23 1WA Groot-Brittannië Tel: +44 (1) 670 734 093 Fax: +44 (1) 670 732 537 E-mail: customerservice.bioprocessing@thermofisher.com

9.3 Technische ondersteuning

Technische ondersteuning voor de HyPerforma S.U.M.'s is beschikbaar in verschillende indelingen. Een gedeelte of alles van het volgende kan van toepassing zijn, afhankelijk van persoonlijke ervaringen en omstandigheden.

Hotline en e-mail voor technisch onderhoud

Neem contact op met uw verkoopvertegenwoordiger van Thermo Fisher Scientific voor algemene productprijzen, beschikbaarheid, levering, bestelinformatie en klachten over producten.

Bel +1 435 792 8500 (Verenigde Staten) of +44 (1) 670 734 093 (Europa, Verenigd Koninkrijk) voor directe en onmiddellijke antwoorden op algemene vragen over producten en technische productinformatie (technische ondersteuning). U kunt ook contact opnemen met de technische ondersteuning door een e-mail te sturen naar: techsupport.bioprocessing@thermofisher.com

Eerste installatie en gebruik

De juiste technische ondersteuning is beschikbaar om te helpen bij de eerste installatie van ieder S.U.M.-systeem. Als u hulp nodig heeft bij het installeren en gebruiken van uw S.U.M.-systeem, vraag hier dan naar als u het product koopt.

Training

Er zijn trainingen beschikbaar voor het opstarten en gebruiken van de S.U.M. Neem voor meer informatie contact op met uw verkoopvertegenwoordiger van Thermo Fisher Scientific.

thermoscientific

Bekijk meer informatie op thermofisher.com/sut

Voor onderzoek of verdere productie. Niet voor diagnostisch gebruik of directe toedieningen in mensen of dieren.

© 2021 Thermo Fisher Scientific Inc. Alle rechten voorbehouden. Alle handelsmerken zijn het eigendom van Thermo Fisher Scientific en haar dochterondernemingen tenzij anders vermeld. CPC en AseptiQuik zijn handelsmerken van Colder Products Company. Pall, Kleenpak en Acro zijn handelsmerken van Pall Corporation. Mettler Toledo is een handelsmerk van Mettler Toledo AG. SmartSite is een handelsmerk van Becton, Dickinson and Company. Teflon is een handelsmerk van The Chemours Company. SterilEnz is een handelsmerk van PAW BioScience. PendoTECH is een handelsmerk van PendoTECH. AC-Tech is een handelsmerk van Lenze. PharMed, PharmaPure en C-Flex zijn handelsmerken van Saint Gobain Performance Plastics. BioQuate is een handelsmerk van General Electric Company. Gore is een handelsmerk van W. L. Gore & Associates, Inc. Lynx is een handelsmerk van EMD Millipore Corporation. Watson-Marlow is een handelsmerk van Watson-Marlow, Inc. Clave is een handelsmerk van Victus. PreSens is een handelsmerk van PreSens Precision Sensing GmbH. **DOC0042NL Rev. C 0321**

