

# Thermo Scientific Solaris -ravistinsarja

# Käyttöohjeet

70900204-е • 2024-11



# Sisältö

Esipuhe		5
Tietoja tä	stä käsikirjasta	5
Kävttötar	koitus	5
,, <b>,</b>	Avoin ravistin	5
	Inkuboitu ja jäähdyttävä ravistin	5
	Tarvittavat taidot	5
Signaalisa	anat ja symbolit	6
Ū	Ravistimessa ja lisävarusteissa käytetyt symbolit	6
	Käyttöohjeessa käytetyt symbolit	7
Turvaohje	et	7
1 Teknico	t tiedet (	
I. Teknise		
1.1. Tek	niset tiedot	11
	Thermo Scientific Solaris 2000 ravistin	11
	Thermo Scientific Solaris 4000 ravistin	12
	Thermo Scientific Solaris 2000 I ravistin	13
	Thermo Scientific Solaris 4000 I ravistin	14
	Thermo Scientific Solaris 2000 R ravistin	15
	Thermo Scientific Solaris 4000 R ravistin	16
1. 2. Tarv	/ikkeet	17
1. 2. 1.	Alustat	18
1. 2. 2.	Koepullon pidikkeet	19
1. 2. 3.	Mikrolevyn / DeepWell-levyn pidikkeet	23
1. 2. 4.	Koeputkitelineet	24
1. 2. 5.	Kallistuskulmaltaan säädettävien koeputkitelineiden	~~
ріаіккеє		28
1.2.6.		30
1. 2. 7.	Nelikulmaisen reagenssipullon pidikkeet	32
1. 2. 8.	Nalgene-dekantterilasin telineet	34
1. 2. 9.		38
1. 2. 10.		39
1. 2. 11.	Liimakiinnitteiset matot	40
1. 2. 12.	. Yleiset lisätarvikkeet	42

1. 3. Direktiivit ja	ı standardit	43
1.4. Tuotekatsau	JS	44
1. 4. 1. Solaris 2	.000	44
1. 4. 2. Solaris 4	.000	44
1. 4. 3. Solaris 2	2000 I / 2000 R	45
1. 4. 4. Solaris 4	.000 I / 4000 R	45
1. 4. 5. Liitännät		46
2. Kuljetus ja asei	nnus	49
2. 1. Purku		49
Tuotteet		49
2. 2. Sijoituspaik	ka	50
2.3. Kulietus		52
Ravistim	ien käsittelv	52
2 4 Vaaitus	······································	53
2.5. Verkkovirtal	liitäntä	53
		55
2. 6. Ensimmaine	an kaynnistys	54
2.7. Varastointi .		59
2.8. Toimitus		59
3. Käyttö		60
3.1. Virta päälle	/ pois	60
3 2 Graafinen k	ävttöliittymä	60
3. 2. 1. Ravistim	en käytön perusparametrien määritys	61
3. 2. 2. Tila		66
3. 2. 3. Asetukse	ət	68
3. 2. 4. Ohjelmat	t	78
3.3. Tarvikkeet .		88
3. 3. 1. Alustan a	asennus	89
3. 3. 2. Koepullo	pidikkeen ja astian asennus	92
3. 3. 3. Nelikulm	aisen pidikkeen asennus	94
3.3.4. Koeputki	itelineen asennus	96
3.3.5. Kallistus	kulmaltaan säädettävän koeputkitelineen	
pidikkeen asennu	IS	97
3. 3. 6. Mikrolev	yn / Deepwell-levyn pidikkeen asennus	98

3. 3. 7. Säädettävän astianpidikkeen asennus	. 99
3. 3. 8. Dekantterilasitelineen asennus	101
3. 3. 9. Erotussuppilon pidikkeen asennus	102
3. 3. 10. Tarviketason asennus	105
3. 3. 11. Liimakiinnitteisen maton asennus ja käyttö	106
3. 3. 12. Kaasunjakoputkiston asennus	113
3. 4. Lastaus ja normaali käyttö	115
Lataa	116
Normaali käyttö	117
4. Huolto ja kunnossapito	121
4.1. Perusteita	122
Apuvälineiden tarkistus	123
4. 2. Puhdistus.	123
Kosketusnäyttö	123
4. 3. Desinfiointi	124
4. 4. Dekontaminaatio	124
4. 5. Autoklavointi	125
4. 6. Lämpötilan kalibrointi	126
4. 6. 1. Lämpötilan kalibroinnin parhaat käytännöt	126
4. 6. 2. Lämpötilan kalibrointimenettely	126
4. 6. 3. Lämpötilan kalibroinnin tarkistus	130
4. 7. Laiteohjelmiston asennus	130
4. 8. Alustan vaihto	134
4.9. Kunnossapito	134
4. 10. Toimitus ja hävittäminen	134
5. Vianmääritys	135
GPL (yleinen julkinen lisenssi)	137
Sisältö	138
Olgano	100

# Esipuhe

# Tietoja tästä käsikirjasta

Tässä käyttöoppaassa kuvataan seuraavat ravistimet ja yhteensopivat lisävarusteet ("1. 2. Tarvikkeet" sivulla 17):

Ravistin	Tuotenro
Thermo Scientific Solaris 2000 ravistin, 100–240 V $\pm 10$ %, 50 / 60 Hz	SK2000
Thermo Scientific Solaris 4000 ravistin, 100–240 V $\pm 10$ %, 50 / 60 Hz	SK4000
Thermo Scientific Solaris 2000 I ravistin, 100–120, 200–240 V $\pm 10~\%,$ 50 / 60 Hz	SK2001
Thermo Scientific Solaris 4000 I ravistin, 100–120, 200–240 V $\pm 10~\%,$ 50 / 60 Hz	SK4001
Thermo Scientific Solaris 2000 R ravistin, 100–240 V $\pm 10$ %, 50 / 60 Hz	SK2002
Thermo Scientific Solaris 4000 R ravistin, 100-240 V ±10 %, 50 / 60 Hz	SK4002

Tämän käyttöoppaan kuvauksissa ja ohjeissa viitataan kuhunkin ravistimeen sen luokan mukaisesti eikä ravistimen täydellisellä tuotenimellä:

- Avoin ravistin
- Inkuboitu ravistin
- Jäähdyttävä ravistin

# Käyttötarkoitus

### Avoin ravistin

Laite, jota käytetään tutkimuksessa ja valmistuksessa ravistamaan liuoksia 2D-radalla samalla kun kontrolloidaan nopeutta ja aikaa sovelluksissa.

### Inkuboitu ja jäähdyttävä ravistin

Laite, jota käytetään tutkimuksessa ja valmistuksessa liuosten ravistukseen 2D-radalla ja joka samalla säätelee nopeutta, aikaa ja lämpötilaa kulloisessakin käyttötilanteessa.

### Tarvittavat taidot

Tätä ravistinta saa käyttää vain koulutettu henkilökunta.

# Signaalisanat ja symbolit

Signaalisana	Vaaran aste
VAROITUS	Osoittaa vaarallisen tilanteen, joka voi johtaa kuolemaan tai vakavaan vammaan, jos sitä ei vältetä.
NOUDATA VAROVAISUUTTA	Näyttää vaarallisen tilanteen, jota välttymättä voi seurata pieni tai keskisuuri vamma.
ниоміо	Näyttää tietoja, jotka arvioidaan tärkeiksi, mutta eivät vaaraa aiheuttaviksi.

# Ravistimessa ja lisävarusteissa käytetyt symbolit

Huomioi ohjekäsikirjassa olevat ohjeet pitääksesi itsesi turvassa ja ympäristösi turvallisena.

Seuraa käyttöopasta
Irrota pääpistoke
Yleinen vaara
Kuuman pinnan aiheuttama vaara
Mekaanisten osien sulkeutuminen vahingossa voi aiheuttaa sormi- tai käsivamman

# Käyttöohjeessa käytetyt symbolit

Huomioi ohjekäsikirjassa olevat ohjeet pitääksesi itsesi turvassa ja ympäristösi turvallisena.

Yleinen vaara	<u>A</u>	Sähköiskun vaara
Biologinen vaara		Viiltovaara
Vaara syttyvistä materiaaleista	i	Näyttää tietoja, jotka arvioidaan tärkeiksi, mutta eivät vaaraa aiheuttaviksi.
Mekaanisten osien sulkeutuminen vahingossa voi aiheuttaa sormi- tai käsivamman.		

# Turvaohjeet



VAROITUS

Näiden turvaohjeiden noudattamatta jättäminen voi johtaa vaarallisiin tilanteisiin, jotka voivat aiheuttaa kuoleman tai vakavan vammautumisen, jos niitä ei vältetä.

- Seuraa turvaohjeita.
  - Ravistinta saa käyttää vain käyttötarkoitukseensa. Väärä käyttö voi johtaa vammautumiseen, kontaminaatioon ja vammoihin, joilla on kohtalokkaat seuraukset.
  - Ravistinta saa käyttää vain koulutettu henkilökunta.
  - On käyttäjän vastuulla varmistaa, että asianmukaisia suojavarusteita käytetään. Seuraa the World Health Organization:in (WHO) "Laboratorion bioturvallisuuskäsikirjaa" ja maakohtaisia määräyksiä.



#### Vahingot väärästä virtalähteestä.

Pidä huoli että ravistin on kytketty vain kunnolla maadoitettuihin pistorasioihin.

VAROITUS



VAROITUS

#### Riski vaarallisten aineiden käsittelystä.

Kun tehdään työtä syövyttävien aineiden kanssa (suolaliuokset, hapot, emäkset), lisävarusteet ja ravistin täytyy puhdistaa perusteellisesti.

- Ravistin ei ole inertti tai suojattu räjähdyksiä vastaan. Älä koskaan käytä ravistinta räjähdysherkässä ympäristössä.
- Älä sekoita myrkyllisiä tai radioaktiivisia materiaaleja tai mitään patogeenisiä mikro-organismeja ilman sopivia turvatoimenpiteitä.
- Mmitä tahansa vaarallisia materiaaleja sekoitettaessa on huomioitava Maailman terveysjärjestön (WHO) "Laboratorion bioturvallisuuskäsikirja" ja paikalliset säädökset. Kun sekoitetaan mikrobiologisia näytteitä riskiryhmässä II (Maailman terveysjärjestön (WHO) "Laboratorion bioturvallisuuskäsikirjan" mukaan), on käytettävä aerosolivarmoja biologisia tiivisteitä. Katso internetistä sivua World Health Organization (www. who.int) löytääksesi "Laboratory Biosafety Manual'in". Korkeamman riskin ryhmien materiaaleille on käytettävä erikoisturvatoimenpiteitä.
- Jos toksiinit tai patogeeniset aineet ovat kontaminoineet ravistimen tai sen osat, tulee suorittaa asianmukaiset desinfiointi- ja dekontaminointitoimenpiteet ("Dekontaminaatio" sivulla 124; "Desinfiointi" sivulla 124).
- Jos vaarallinen tilanne sattuu, käännä virta pois ravistimesta ja poistu välittömästi alueelta.



VAROITUS

Terveysvahingot tartuntavaarallisista aineista.

Jos nesteitä tai muita aineita vahingossa läikkyy alustan alle, sammuta ravistin välittömästi, irrota se pistorasiasta ja poista alusta ("Alustat" sivulla 18).

Puhdista läikkynyt materiaali seuraten tavanomaisia laboratoriotoimenpiteitä. Käytä kunnollisia henkilökohtaisia suojaimia.



Terveydelle vaarallinen räjähtävien tai tulenarkojen materiaalien tai aineiden ravistelu.

Älä sekoita räjähtäviä tai tulenarkoja materiaaleja tai aineita.

VAROITUS



NOUDATA VAROVAISUUTTA

#### Kuumien pintojen aiheuttamat palovammat.

Lämpötilasäädellyssä ravistimessa pinnat, kuten alusta ja lisävarusteet, voivat kuumentua kannen alla. Älä koske kuumiin pintoihin. Odota, kunnes kuumat pinnat ovat jäähtyneet.



VAROVAISUUTTA

# NOUDATA

#### Viiltohaavat lasinsirpaleista.

Irrallaan oleva alusta ja lisävarusteet, kuten puristimet, voivat johtaa rikkinäiseen lasiin jos astiat putoavat ravistimesta. Varmista että alusta ja lisävarusteet on asianmukaisesti asennettu käyttäen oikeita työkaluja ja ruuveja. Varmista, että puristimet ovat riittävät kantamaan astian ja näytteen painon valitulla nopeudella. Puristimet, joilla on yksilöllisesti konfiguroitu kuorma, voivat omata alhaisemman nopeuskapasiteetin kuin ravistimen stabiliteettirajoitus. Katso "3. 4. Lastaus janormaali käyttö" sivulla 115.

Huomioi oudot äänet. Tämä voi johtua irronneesta alustasta tai lisävarusteesta.



# Ravistimen vaurio tai vaurioituneen kosketusnäytön aiheuttama toimintahäiriö.

NOUDATA VAROVAISUUTTA

- Alä käytä ravistinta, jos kosketusnäyttö on vaurioitunut.
- Katkaise ravistimen virta. Irrota päävirtajohto. Vaihdatuta kosketusnäyttö valtuutetulla huoltomiehellä.



# Biologinen vahinko johtuen rikkinäisistä tai vuotavista astioista.

NOUDATA VAROVAISUUTTA Väärin asennetut apuvälineet voivat johtaa läikkyneisiin näytteisiin.

- Varmista, että apuvälineet on oikein asennettu käyttäen oikeita työkaluja ja ruuveja.
- Pidä huolta, että apuvälineet sopivat alustaan mielekkäästi.
- Käytä aina sopivan kokoista astiaa ja lisävarustetta.
- Astioiden täytyy olla virheettömiä ja oikein asennettu.



# Turvallisuus voi vaarantua väärästä lastaamisesta ja rikkonaisista apuvälineistä.

- NOUDATA VAROVAISUUTTA
- Varmista aina että kuorma (apuvälineet ja näytteet) on jakautunut niin tasaisesti kuin mahdollista, erityisesti kun käytät kaksoispinoalustaa.
- Älä käytä apuvälineitä, joissa on merkkejä korroosiosta tai halkeamia. Ota yhteys asiakaspalveluun lisätietoja varten.
- Käytä vain ravistimia, jotka on oikein lastattu.
- Älä ylikuormita ravistinta.
- Varmista, että kaikki apuvälineet on asennettu oikein ennen ravistimen käyttöä. Noudata ohjeita osiossa "Tarvikkeet" sivulla 88.



# Perusasioiden jättäminen huomiotta ja sen tuoma fyysinen vahinko.

- Älä käytä ravistinta ilman kunnolla asennettua alustaa.
- Älä käytä ravistinta, jos siitä puuttuu osia tai sen ulkopuoli on vaurioitunut tai se puuttuu.
- Älä käytä ravistinta, jonka kannen jousi on viallinen.
  Viallinen kannen jousi ei pysty pitämään kantta turvallisesti täysin auki.
- Älä siirrä ravistinta kun se käy.
- Älä nojaa ravistimeen.
- Älä kuormita tai pura ravistinta ennen kuin se on pysähtynyt täysin ja tämä on vahvistettu kosketusnäytöllä.
- Älä pane mitään ravistimen päälle sen ollessa käynnissä.
- Älä kosketa alustaa tai mitään apuvälineitä ravistimessa sen ollessa käynnissä.
- Ravistimen koteloa ei saa avata käyttäjän toimesta.



# Suojaus saattaa heikentyä sopimattomien apuvälineiden takia.

# ниоміо

Käytä ravistimessa vain Thermo Fisher Scientificin hyväksymiä apuvälineitä. Päivitettyjä listoja varten tarkista www.thermofisher.com.



HUOMIO

### Sammuttaaksesi ravistimen:

Paina the STOP-kytkintä. Käännä ravistin pois päältä pääkytkimestä. Irrota sähkökosketin seinästä. Hätätilassa kytke virta pois virtakytkimestä.

#### NOUDATA VAROVAISUUTTA

# 1. Tekniset tiedot

# 1.1. Tekniset tiedot

Thermo Scientific Solaris 2000 ravistin				
Nopeusvalikoima	15–525 rpm			
Ajoaika	99 h 59 min (1 min lisäyksellä) tai jatkuva moodi			
Melutaso maksiminopeudella	50 dB (A) (1 m laitteen e	edessä 1,6 m korkeudella)		
Maksimikuorma (sis. alusta, apuvälineet ja näytteet)	25 kg (55 lbs)	25 kg (55 lbs)		
Sähkökytkentä	100–240 V ±10 %, 50 /	′ 60 Hz		
Virrankulutus	60 W			
Ympäristöolosuhteet				
Säilytystä ja kuljetusta varten	Lämpötila: -10 °C - 55 ° Kosteus: 15 % 85 %	°C		
Käyttöä varten	Käyttö sisätiloissa Korkeudet 3000:een metriin saakka merenpinnan yläpuolella. Lämpötila: 5 °C - 40 °C Maks. suhteellinen kosteus 80 %, 31 °C lämpötilaan asti.; laskee lineaarisesti suhteelliseen kosteuteen 50 %; 40 °C lämpötilassa.			
Saastuttamisaste	2			
Ylijännitekategoria	1			
IP	20			
Liittymät	USB	Ethernet		
	2x USB-A 2.0	RJ45		
Mitat	Pituus	Leveys	Korkeus	
	47 cm (18,5 in)	37 cm (14,5 in)	15 cm (5,5 in)	
Paino	20,9 kg (46,0 lbs)			

Taulukko 1: Tekniset tiedot Solaris 2000

Nopeusvalikoima	15–525 rpm		
Ajoaika	99 h 59 min (1 min lisäyksellä) tai jatkuva moodi		
Melutaso maksiminopeudella	54 dB (A) (1 m laitteen e	dessä 1,6 m korkeudella)	
Maksimikuorma (sis. alusta, apuvälineet ja näytteet)	43 kg (95 lbs)		
Sähkökytkentä	100–240 V ±10 %, 50 /	60 Hz	
Virrankulutus	80 W		
Ympäristöolosuhteet			
Säilytystä ja kuljetusta varten	Lämpötila: -10 °C - 55 ° Kosteus: 15 % 85 %	С	
Käyttöä varten	Käyttö sisätiloissa Korkeudet 3000:een metriin saakka merenpinnan yläpuolella. Lämpötila: 5 °C - 40 °C Maks. suhteellinen kosteus 80 %, 31 °C lämpötilaan asti.; laskee lineaarisesti suhteelliseen kosteuteen 50 %; 40 °C lämpötilassa.		
Saastuttamisaste	2		
Ylijännitekategoria	1		
P	20		
Liittymät	USB	Ethernet	
	2x USB-A 2.0	RJ45	
Mitat	Pituus	Leveys	Korkeus
	65 cm (25,6 in)	58 cm (22,8 in)	18 cm (7 in)
Paino	75,1 kg (165,5 lbs)		

Taulukko 2: Tekniset tiedot Solaris 4000

# Thermo Scientific Solaris 2000 I ravistin

Nopeusvalikoima	15–525 rpm		
Lämpötila-asetuksen arvorajat	30-60 °C		
Koepullon lämpötilan stabiilius lämpötilassa 37 °C	Ympäristön lämpötila 23 °C; 1 tunnin ajan, kun kammion lämpötila on vakaa $\pm$ 0,1 °C		
Lämpötilan tasaisuus koepullossa lämpötilassa 37 °C	Ympäristön lämpötila 23 ° ± 0,5 °C	'C; 1 tunnin ajan, kun kamn	nion lämpötila on vakaa
Ajoaika	99 h 59 min (1 min lisäyl	ksellä) tai jatkuva moodi	
Melutaso maksiminopeudella	52 dB (A) (1 m laitteen e	dessä 1,6 m korkeudella)	
Maksimikuorma (sis. alusta, apuvälineet ja näytteet)	16 kg (35 lbs), mukaan lukien kokoa 11	x 14 oleva alusta painolla	2,3 kg (5,1 lbs)
Sähkökytkentä	100–120, 200–240 V ±	10 %, 50 / 60 Hz	
Virrankulutus	900 W		
Ympäristöolosuhteet			
Säilytystä ja kuljetusta varten	Lämpötila: -10 °C - 55 °C Kosteus: 15 % 85 %		
Käyttöä varten	Käyttö sisätiloissa Korkeudet 3000:een metriin saakka merenpinnan yläpuolella. Lämpötila: 5 °C - 40 °C Maks. suhteellinen kosteus 80 %, 31 °C lämpötilaan asti.; laskee lineaarisesti suhteelliseen kosteuteen 50 %; 40 °C lämpötilassa.		
Saastuttamisaste	2		
Ylijännitekategoria			
P	20		
Liittymät	USB	Ethernet	
	2x USB-A 2.0	RJ45	
Mitat	Pituus	Leveys	Korkeus
	70 cm (27,6 in)	36 cm (14,2 in)	46 cm (18,1 in)
Kansi auki			79 cm (30,8 in)
Paino	44,9 kg (99 lbs)		

Taulukko 3: Tekniset tiedot Solaris 2000 I

#### Tekniset tiedot

# Thermo Scientific Solaris 4000 I ravistin

Nopeusvalikoima	15–525 rpm		
Lämpötila-asetuksen arvorajat	30–60 °C		
Koepullon lämpötilan stabiilius lämpötilassa 37 °C	Ympäristön lämpötila 23 °C; 1 tunnin ajan, kun kammion lämpötila on vakaa $\pm$ 0,1 °C		
Lämpötilan tasaisuus koepullossa lämpötilassa 37 °C	Ympäristön lämpötila 23 ° ± 0,5 °C	°C; 1 tunnin ajan, kun kamn	nion lämpötila on vakaa
Ajoaika	99 h 59 min (1 min lisäyl	ksellä) tai jatkuva moodi	
Melutaso maksiminopeudella	52 dB (A) (1 m laitteen e	dessä 1,6 m korkeudella)	
Maksimikuorma (sis. alusta, apuvälineet ja näytteet)	32 kg (70,5 lbs), mukaan lukien kokoa 18	x 18 oleva alusta painolla	4,3 kg (9,5 lbs)
Sähkökytkentä	100–120, 200–240 V ±	10 %, 50 / 60 Hz	
Virrankulutus	900 W		
Ympäristöolosuhteet			
Säilytystä ja kuljetusta varten	Lämpötila: -10 °C - 55 °C Kosteus: 15 % 85 %		
Käyttöä varten	Käyttö sisätiloissa Korkeudet 3000:een metriin saakka merenpinnan yläpuolella. Lämpötila: 5 °C - 40 °C Maks. suhteeliinen kosteus 80 %, 31 °C lämpötilaan asti.; laskee lineaarisesti suhteeliiseen kosteuteen 50 %; 40 °C lämpötilassa.		
Saastuttamisaste	2		
Ylijännitekategoria			
P	20		
Liittymät	USB	Ethernet	
	2x USB-A 2.0	RJ45	
Mitat	Pituus	Leveys	Korkeus
	77 cm (30,3 in)	57 cm (22,4 in)	55 cm (21,7 in)
Kansi auki			94 cm (36,7 in)
Paino	68,1 kg (150,1 lbs)		

Taulukko 4: Tekniset tiedot Solaris 4000 I

# Thermo Scientific Solaris 2000 R ravistin

Nopeusvalikoima	15–525 rpm			
Lämpötila-asetuksen arvorajat	5–60 °C			
Koepullon lämpötilan stabiilius lämpötilassa 37 °C	Ympäristön lämpötila 23 °C; 1 tunnin ajan, kun kammion lämpötila on vakaa $\pm$ 0,1 °C			
Lämpötilan tasaisuus koepullossa lämpötilassa 37 °C	Ympäristön lämpötila 23 ° ± 0,5 °C	'C; 1 tunnin ajan, kun kamm	nion lämpötila on vakaa	
Ajoaika	99 h 59 min (1 min lisäył	ksellä) tai jatkuva moodi		
Melutaso maksiminopeudella	56 dB (A) (1 m laitteen e	dessä 1,6 m korkeudella)		
Maksimikuorma (sis. alusta, apuvälineet ja näytteet)	16 kg (35 lbs), mukaan lukien kokoa 11	x 14 oleva alusta painolla	2,3 kg (5,1 lbs)	
Sähkökytkentä	100–240 V ±10 %, 50 /	60 Hz		
Virrankulutus	350 W			
Ympäristöolosuhteet				
Säilytystä ja kuljetusta varten	Lämpötila: -10 °C - 55 °C Kosteus: 15 % 85 %			
Käyttöä varten	Käyttö sisätiloissa Korkeudet 3000:een metriin saakka merenpinnan yläpuolella. Lämpötila: 5 °C - 40 °C Maks. suhteellinen kosteus 80 %, 31 °C lämpötilaan asti.; laskee lineaarisesti suhteelliseen kosteuteen 50 %; 40 °C lämpötilassa.			
Saastuttamisaste	2			
Ylijännitekategoria	ll			
P	20			
Liittymät	USB	Ethernet		
	2x USB-A 2.0	RJ45		
Mitat	Pituus	Leveys	Korkeus	
	70 cm (27,6 in)	36 cm (14,2 in)	46 cm (18,1 in)	
Kansi auki			79 cm (30,8 in)	
Paino	47,2 kg (104,1 lbs)			

Taulukko 5: Tekniset tiedot Solaris 2000 R

# Thermo Scientific Solaris 4000 R ravistin

Nopeusvalikoima	15–525 rpm				
Lämpötila-asetuksen arvorajat	4–60 °C	4–60 °C			
Koepullon lämpötilan stabiilius lämpötilassa 37 °C	Ympäristön lämpötila 23 °C; 1 tunnin ajan, kun kammion lämpötila on vakaa $\pm$ 0,1 °C				
Lämpötilan tasaisuus koepullossa lämpötilassa 37 °C	Ympäristön lämpötila 23 °C ± 0,5 °C	C; 1 tunnin ajan, kun kammi	on lämpötila on vakaa		
Ajoaika	99 h 59 min (1 min lisäyks	sellä) tai jatkuva moodi			
Melutaso maksiminopeudella	56 dB (A) (1 m laitteen ed	essä 1,6 m korkeudella)			
Maksimikuorma (sis. alusta, apuvälineet ja näytteet)	32 kg (70,5 lbs), mukaan lukien kokoa 18 >	< 18 oleva alusta painolla 4	,3 kg (9,5 lbs)		
Sähkökytkentä	100–240 V ±10 %, 50 / 6	60 Hz			
Virrankulutus	750 W				
Ympäristöolosuhteet					
Säilytystä ja kuljetusta varten	Lämpötila: -10 °C - 55 °C Kosteus: 15 % 85 %				
Käyttöä varten	Käyttö sisätiloissa Korkeudet 3000:een metriin saakka merenpinnan yläpuolella. Lämpötila: 5 °C - 40 °C Maks. suhteellinen kosteus 80 %, 31 °C lämpötilaan asti.; laskee lineaarisesti suhteelliseen kosteuteen 50 %; 40 °C lämpötilassa.				
Saastuttamisaste	2				
Ylijännitekategoria	11				
IP	20				
Liittymät	USB	Ethernet			
	2x USB-A 2.0	RJ45			
Mitat	Pituus	Leveys	Korkeus		
	77 cm (30,3 in)	57 cm (22,4 in)	55 cm (21,7 in)		
Kansi auki			94 cm (36,7 in)		
Paino	74,9 kg (165,1 lbs)				

Taulukko 6: Tekniset tiedot Solaris 4000 R

# 1.2. Tarvikkeet



Suojaus saattaa heikentyä sopimattomien apuvälineiden takia.

NOUDATA VAROVAI-SUUTTA Käytä ravistimessa vain Thermo Fisher Scientificin hyväksymiä apuvälineitä.

Päivitettyjä listoja varten tarkista www.thermofisher.com.

# 1.2.1. Alustat

Alusta		
Solaris 2000	Paino	
Thermo Scientific Solaris 12x14 yleisalusta	2,7 kg (5,9 lbs)	SK1214
Thermo Scientific Solaris 12x14 Dual Stack yleisalusta	6,4 kg (14,1 lbs)	SK1214D
Thermo Scientific Solaris 12x14 Dual Stack Yleisalusta päivityssarja		SK1214DK
Thermo Scientific Solaris 18x18 yleisalusta	4,3 kg (9,4 lbs)	SK1818
Thermo Scientific Solaris 18x18 Dual Stack yleisalusta	9,7 kg (21,3 lbs)	SK1818D
Thermo Scientific Solaris 18x18 Dual Stack Yleisalusta päivityssarja		SK1818DK
Thermo Scientific Solaris 18x24 yleisalusta	6,0 kg (13,2 lbs)	SK1824
Solaris 4000		
Thermo Scientific Solaris 18x30 yleisalusta	7,5 kg (16,5 lbs)	SK1830
Thermo Scientific Solaris 18x30 Dual Stack yleisalusta	16,3 kg (35,9 lbs)	SK1830D
Thermo Scientific Solaris 18x30 Dual Stack Yleisalusta päivityssarja		SK1830DK
Thermo Scientific Solaris 36x24 yleisalusta	12,0 kg (26,4 lbs)	SK3624
Solaris 2000 I / 2000 R		
Thermo Scientific Solaris 11x14 yleisalusta	2,3 kg (5,1 lbs)	SK1114
Solaris 4000 I / 4000 R		
Thermo Scientific Solaris 18x18 yleisalusta	4,3 kg (9,5 lbs)	SK1818
Varaosasarjat ja lisävarusteet		
Puristimen varaosasarja (Ruuveja)		SK1001
Alustan varaosasarja, pieni (alustan ruuvit, työkalu)		SK0100
Alustan varaosasarja, suuri (alustan ruuvit, työkalu)		SK0101
Ruuvitaltta lisävarusteille		75004131

Taulukko 7: Saatavissa olevat alustat

# 1. 2. 2. Koepullon pidikkeet



## Solaris 2000 -alustoille

	Tuotenro		Yleinen		Dual	Stack
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
10 ml Erlenmeyer	30150BI	72	113	157	140	226
25 ml Erlenmeyer	30151	42	64	80	80	124
50 ml Erlenmeyer	30152BI	42	64	80	80	124
125 ml Erlenmeyer	30153	15	32	40	30	62
250 ml Erlenmeyer	30154BI	9	16	24	16	32
300 ml Erlenmeyer	30155	9	16	20	16	32
500 ml Erlenmeyer	30156BI	9	16	20	16	32
1   Erlenmeyer	30157BI	4	9	11	8	16
2 l Erlenmeyer	30158	3	5	6	-	-
4 l Erlenmeyer	30159	1	4	4	-	-
5 l Erlenmeyer	30159B	1	2	4	-	-
6 l Erlenmeyer	30160	1	2	2	-	-
2 800 ml Fernbach Flask	30162	1	4	4	-	-
Low form viljelmäpullo 2,5 l	30161	1	2	2	-	-

Taulukko 8: Saatavilla olevat koepullon pidikkeet Solaris 2000 -alustoille

# Solaris 4000 -alustoille

	Tuotenro	Yle	inen	Dual Stack
		18 x 30	36 x 24	18 x 30
10 ml Erlenmeyer	30150BI	203	187	402
25 ml Erlenmeyer	30151	112	187	220
50 ml Erlenmeyer	30152BI	112	187	220
125 ml Erlenmeyer	30153	46	83	92
250 ml Erlenmeyer	30154BI	34	40	68
300 ml Erlenmeyer	30155	28	40	56
500 ml Erlenmeyer	30156BI	28	40	56
1   Erlenmeyer	30157BI	14	20	28
2 l Erlenmeyer	30158	8	11	16
4 l Erlenmeyer	30159	6	8	-
5 l Erlenmeyer	30159B	5	8	-
6 l Erlenmeyer	30160	3	6	-
2 800 ml Fernbach Flask	30162	6	8	12
Low Form viljelmäpullo 2,5 I	30161	3	6	6

Taulukko 9: Saatavilla olevat koepullon pidikkeet Solaris 4000 -alustoille

## Solaris 2000 I / 2000 R -alustoille

	Tuotenro	Yleinen
		11 x 14
10 ml Erlenmeyer	30150BI	59
25 ml Erlenmeyer	30151	35
50 ml Erlenmeyer	30152BI	35
125 ml Erlenmeyer	30153	15
250 ml Erlenmeyer	30154BI	7
300 ml Erlenmeyer	30155	7
500 ml Erlenmeyer	30156BI	7
1   Erlenmeyer	30157BI	4
2 l Erlenmeyer	30158	2
4 l Erlenmeyer	30159	-
5 l Erlenmeyer	30159B	-
6 l Erlenmeyer	30160	-
2 800 ml Fembach Flask	30162	1
Low Form viljelmäpullo 2,5 l	30161	1

Taulukko 10: Saatavilla olevat koepullon pidikkeet Solaris 2000 I / 2000 R -alustalle

## Solaris 4000 I / 4000 R -alustoille

	Tuotenro	Yleinen
		18 x 18
10 ml Erlenmeyer	30150BI	113
25 ml Erlenmeyer	30151	64
50 ml Erlenmeyer	30152BI	64
125 ml Erlenmeyer	30153	32
250 ml Erlenmeyer	30154BI	16
300 ml Erlenmeyer	30155	16
500 ml Erlenmeyer	30156BI	16
1   Erlenmeyer	30157BI	9
2 l Erlenmeyer	30158	6
4 l Erlenmeyer	30159	-
5 l Erlenmeyer	30159B	-
6 l Erlenmeyer	30160	-
2 800 ml Fembach Flask	30162	4
Low Form viljelmäpullo 2,5 l	30161	1

Taulukko 11: Saatavilla olevat koepullon pidikkeet Solaris 4000 I / 4000 R -alustalle

# 1. 2. 3. Mikrolevyn / DeepWell-levyn pidikkeet



#### Solaris 2000 -alustoille

	Tuotenro	Yleinen		Dual Stack		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Mikrolevy / DeepWell-levy	30175	7	12	16	14	24

Taulukko 12: Mikrolevyn/DeepWell-levyn pidikkeiden kapasiteetti Solaris 2000 -alustoilla

# Solaris 2000 I / 2000 R -alustoille

	Tuotenro	Yleinen
		11 x 14
Mikrolevy / DeepWell-levy	30175	6

Taulukko 13: Mikrolevyn/DeepWell-levyn pidikkeiden kapasiteetti Solaris 2000 I / 2000 R -alustalla

### Solaris 4000 -alustoille

	Tuotenro	Ylei	Yleinen	
		18 x 30	36 x 24	18 x 30
Mikrolevy / DeepWell-levy	30175	23	36	42

Taulukko 14: Mikrolevyn/DeepWell-levyn pidikkeiden kapasiteetti Solaris 4000 -alustalla

### Solaris 4000 I / 4000 R -alustoille

	Tuotenro	Yleinen
		18 x 18
Mikrolevy / DeepWell-levy	30175	10

Taulukko 15: Mikrolevyn/DeepWell-levyn pidikkeiden kapasiteetti Solaris 4000 I / 4000 R -alustalla

# 1. 2. 4. Koeputkitelineet

#### Solaris 2000 -alustoille



	Tuotenro		Yleinen		Dual	Stack
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Puolikas koko						
10–13 mm, punainen, 6 x 6 teline	30181	8	12	15	14	22
14–16 mm, oranssi, 6 x 6 teline	30183	5	9	11	9	16
17–20 mm, valkoinen, 4 x 5 teline	30185	7	11	14	12	20
21–25 mm, sininen, 4 x 4 teline	30187	6	9	11	10	16
26–30 mm, vihreä, 3 x 3 teline	30189	6	9	12	11	17
Mikrosentrifugi, 1,5 ml, sininen, 4 x 6 teline	30191	6	10	13	11	19
Täysi koko						
10–13 mm, 6 x 12 teline	30180BI	3	7	9	6	12
14–16 mm, 6 x 12 teline	30182	3	4	6	6	8
17–20 mm, 4 x 10 teline	30184	3	5	7	6	8
21–25 mm, 4 x 10 teline	30186	2	3	5	4	6
26–30 mm, 3 x 8 teline	30188	3	4	5	5	8
1,5 ml mikrosentrifugi, 8 x 12 teline	30190	3	4	7	6	8

Taulukko 16: Saatavana olevat koeputkitelineet Solaris 2000-alustoille

# Solaris 4000 -alustoille

	Tuotenro	Yleinen		Dual Stack	
		18 x 30	36 x 24	18 x 30	
Puolikas koko					
10–13 mm, punainen, 6 x 6 teline	30181	21	32	40	
14–16 mm, oranssi, 6 x 6 teline	30183	13	20	24	
17–20 mm, valkoinen, 4 x 5 teline	30185	18	20	34	
21–25 mm, sininen, 4 x 4 teline	30187	14	22	26	
26–30 mm, vihreä, 3 x 3 teline	30189	15	24	30	
1,5 ml Mikrosentrifugi, sininen, 4 x 6 teline	30191	18	24	34	
Täysi koko					
10–13 mm, 6 x 12 teline	30180BI	13	20	18	
14–16 mm, 6 x 12 teline	30182	9	12	16	
17–20 mm, 4 x 10 teline	30184	9	15	18	
21–25 mm, 4 x 10 teline	30186	6	9	12	
26–30 mm, 3 x 8 teline	30188	7	10	13	
1,5 ml mikrosentrifugi, 8 x 12 teline	30190	9	12	17	

Taulukko 17: Saatavana olevat koeputkitelineet Solaris 4000-alustoille

# Solaris 2000 I / 2000 R -alustoille

	Tuotenro	Yleinen
Puolikas koko		
10–13 mm, punainen, 6 x 6 teline	30181	6
14–16 mm, oranssi, 6 x 6 teline	30183	2
17–20 mm, valkoinen, 4 x 5 teline	30185	4
21–25 mm, sininen, 4 x 4 teline	30187	2
26–30 mm, vihreä, 3 x 3 teline	30189	4
Mikrosentrifugi, 1,5 ml, sininen, 4 x 6 teline	30191	3
Täysi koko		
10–13 mm, 6 x 12 teline	30180BI	3
14–16 mm, 6 x 12 teline	30182	2
17–20 mm, 4 x 10 teline	30184	2
21–25 mm, 4 x 10 teline	30186	1
26–30 mm, 3 x 8 teline	30188	2
1,5 ml mikrosentrifugi, 8 x 12 teline	30190	2

Taulukko 18: Saatavana olevat koeputkitelineet Solaris 2000 I / 2000 R-alustoille

# Solaris 4000 I / 4000 R -alustoille

	Tuotenro	Yleinen
		18 x 18
Puolikas koko		
10–13 mm, punainen, 6 x 6 teline	30181	12
14–16 mm, oranssi, 6 x 6 teline	30183	8
17–20 mm, valkoinen, 4 x 5 teline	30185	10
21–25 mm, sininen, 4 x 4 teline	30187	8
26–30 mm, vihreä, 3 x 3 teline	30189	9
Mikrosentrifugi, 1,5 ml, sininen, 4 x 6 teline	30191	8
Täysi koko		
10–13 mm, 6 x 12 teline	30180BI	7
14–16 mm, 6 x 12 teline	30182	4
17–20 mm, 4 x 10 teline	30184	5
21–25 mm, 4 x 10 teline	30186	3
26–30 mm, 3 x 8 teline	30188	4
1,5 ml mikrosentrifugi, 8 x 12 teline	30190	5

Taulukko 19: Saatavana olevat koeputkitelineet Solaris 4000 I / 4000 R-alustoille

# 1. 2. 5. Kallistuskulmaltaan säädettävien koeputkitelineiden pidikkeet



### Solaris 2000 -alustoille

	Tuotenro	Yleinen			Dual Stack		
	laotonio				D'aa.	otaon	
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18	
10–13 mm, 72 paikkaa	236090	3	4	6	6	8	
16–20 mm, 40 paikkaa	236091	3	4	6	5	8	
21–25 mm, 40 paikkaa	236092	3	3	4	5	6	
26–30 mm, 24 paikkaa	236093	3	3	4	5	6	

Taulukko 20: Saatavilla olevat kallistuskulmaltaan säädettävät koeputkitelinepidikkeet Solaris 2000 -alustoille

### Solaris 4000 -alustoille

	Tuotenro	Yleinen		Dual Stack	
		18 x 30	36 x 24	18 x 30	
10–13 mm, 72 paikkaa	236090	7	14	14	
16–20 mm, 40 paikkaa	236091	7	14	14	
21–25 mm, 40 paikkaa	236092	5	9	10	
26–30 mm, 24 paikkaa	236093	6	9	12	

Taulukko 21: Saatavilla olevat kallistuskulmaltaan säädettävät koeputkitelinepidikkeet Solaris 4000 -alustoille

# Solaris 2000 I / 2000 R -alustoille

	Tuotenro	Yleinen
		11 x 14
10–13 mm, 72 paikkaa	236090	2
16–20 mm, 40 paikkaa	236091	2
21–25 mm, 40 paikkaa	236092	1
26–30 mm, 24 paikkaa	236093	1

Taulukko 22: Saatavilla olevat kallistuskulmaltaan säädettävät koeputkitelinepidikkeet Solaris 2000 I / 2000 R-alustoille

### Solaris 4000 I / 4000 R -alustoille

	Tuotenro	Yleinen
		18 x 18
10–13 mm, 72 paikkaa	236090	4
16–20 mm, 40 paikkaa	236091	3
21–25 mm, 40 paikkaa	236092	3
26–30 mm, 24 paikkaa	236093	3

Taulukko 23: Saatavilla olevat kallistuskulmaltaan säädettävät koeputkitelinepidikkeet Solaris 4000 I / 4000 R-alustoille

# 1. 2. 6. Säädettävät säiliön kiinnikkeet



#### Solaris 2000 -alustoille

	Tuotenro	Yleinen			Dual Stack		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18	
Yhdelle korkeudelle säädettävä astian pidike, 11" x 14"	75004104	-	-	-	-	_	
Kahdelle korkeudelle säädettävä astian pidike, 12" x 14"	75004102	1	1	1	-	2	
Kahdelle korkeudelle säädettävä astian pidike, 18" x 18"	75004103	-	1	1	-	-	
Yhden astian säädettävä pidike	75004101	4	4	6	6	8	

Taulukko 24: Saatavilla olevat astian pidikkeet Solaris 2000 -alustoille

### Solaris 4000 -alustoille

	Tuotenro	Yleinen		Dual Stack	
		18 x 30	36 x 24	18 x 30	
Yhdelle korkeudelle säädettävä astian pidike, 11" x 14"	75004104	-	-	-	
Kahdelle korkeudelle säädettävä astian pidike, 12" x 14"	75004102	2	2	4	
Kahdelle korkeudelle säädettävä astian pidike, 18" x 18"	75004103	1	2	2	
Yhden astian säädettävä pidike	75004101	8	12	16	

Taulukko 25: Saatavilla olevat astian pidikkeet Solaris 4000 -alustoille

# Solaris 2000 I / 2000 R -alustoille

	Tuotenro	Yleinen
		11 x 14
Yhdelle korkeudelle säädettävä astian pidike, 11" x 14"	75004104	1
Kahdelle korkeudelle säädettävä astian pidike, 12" x 14"	75004102	-
Kahdelle korkeudelle säädettävä astian pidike, 18" x 18"	75004103	-
Yhden astian säädettävä pidike	75004101	2

Taulukko 26: Säädettävät astian pidikkeet Solaris 2000 I / 2000 R -alustoille

## Solaris 4000 I / 4000 R -alustoille

	Tuotenro	Yleinen
		18 x 18
Yhdelle korkeudelle säädettävä astian pidike, 11" x 14"	75004104	-
Kahdelle korkeudelle säädettävä astian pidike, 12" x 14"	75004102	1
Kahdelle korkeudelle säädettävä astian pidike, 18" x 18"	75004103	1
Yhden astian säädettävä pidike	75004101	4

Taulukko 27: Säädettävät astian pidikkeet Solaris 4000 I / 4000 R -alustoille

# 1. 2. 7. Nelikulmaisen reagenssipullon pidikkeet



#### Solaris 2000 -alustoille

	Tuotenro	Yleinen		Dual Stack		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
125 ml:n nelikulmaisen pullon pidike	75004106	12	16	24	21	32
250 ml:n nelikulmaisen pullon pidike	75004107	9	16	22	18	32
500 ml:n nelikulmaisen pullon pidike	75004108	6	9	12	11	18
1 000 ml:n nelikulmaisen pullon pidike	75004109	4	9	12	8	18

Taulukko 28: Saatavilla olevat nelikulmaisen reagenssipullon pidikkeet Solaris 2000 -alustoille

#### Solaris 4000 -alustoille

	Tuotenro	Yleinen		Dual Stack
		18 x 30	36 x 24	18 x 30
125 ml:n nelikulmaisen pullon pidike	75004106	28	54	56
250 ml:n nelikulmaisen pullon pidike	75004107	28	52	56
500 ml:n nelikulmaisen pullon pidike	75004108	15	24	30
1 000 ml:n nelikulmaisen pullon pidike	75004109	15	24	28

Taulukko 29: Saatavilla olevat nelikulmaisen reagenssipullon pidikkeet Solaris 4000 -alustoille

# Solaris 2000 I / 2000 R -alustoille

	Tuotenro	Yleinen
		11 x 14
125 ml:n nelikulmaisen pullon pidike	75004106	9
250 ml:n nelikulmaisen pullon pidike	75004107	7
500 ml:n nelikulmaisen pullon pidike	75004108	3
1 000 ml:n nelikulmaisen pullon pidike	75004109	2

Taulukko 30: Saatavilla nelikulmaisen reagenssipullon pidikkeet Solaris 2000 I / 2000 R -alustoille

# Solaris 4000 I / 4000 R -alustoille

	Tuotenro	Yleinen
		18 x 18
125 ml:n nelikulmaisen pullon pidike	75004106	16
250 ml:n nelikulmaisen pullon pidike	75004107	16
500 ml:n nelikulmaisen pullon pidike	75004108	9
1 000 ml:n nelikulmaisen pullon pidike	75004109	9

Taulukko 31: Saatavilla nelikulmaisen reagenssipullon pidikkeet Solaris 4000 I / 4000 R -alustoille

# 1. 2. 8. Nalgene-dekantterilasin telineet

#### Solaris 2000 -alustoille



	Tuotenro		Yleinen		Dual S	Stack
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Koko 11 x 14 / 12 x 14			Те	lineet/paikat	:	
30 ml:n Nalgene-dekantterilasi, 36 paikkaa	75004116	1 / 36	-	-	2/72	-
50 ml:n Nalgene-dekantterilasi, 28 paikkaa	75004129	1 / 28	-	-	2/56	-
100 ml:n Nalgene-dekantterilasi, 16 paikkaa	75004118	1 / 16	-	-	2/32	-
250 ml:n Nalgene-dekantterilasi, 9 paikkaa	75004119	1/9	-	-	2/18	-
400 ml:n Nalgene-dekantterilasi, 8 paikkaa	75004120	1/8	-	-	2/16	-
600 ml:n Nalgene-dekantterilasi, 5 paikkaa	75004121	1/5	-	-	2/10	-
Koko 9 x 18			Те	lineet/paikat	t	
30 ml:n Nalgene-dekantterilasi, 32 paikkaa	75004110	-	2/64	2 / 64	-	4 / 128
50 ml:n Nalgene-dekantterilasi, 24 paikkaa	75004128	-	2 / 48	2/48	-	4 / 96
100 ml:n Nalgene-dekantterilasi, 15 paikkaa	75004112	-	2/30	2/30	-	4 / 60
250 ml:n Nalgene-dekantterilasi, 8 paikkaa	75004113	-	2/16	2/16	-	4 / 32
400 ml:n Nalgene-dekantterilasi, 6 paikkaa	75004114	-	2/12	2/12	-	4 / 24
600 ml:n Nalgene-dekantterilasi, 4 paikkaa	75004115	-	2/8	2/8	-	4 / 16

Taulukko 32: Saatavilla olevat Nalgene-dekantterilasin telineet Solaris 2000 -alustoille

# Solaris 4000 -alustoille

	Tuotenro	Yleinen		Dual Stack
		18 x 30	36 x 24	18 x 30
Koko 11 x 14 / 12 x 14			Telineet/paikat	:
30 ml:n Nalgene-dekantterilasi, 36 paikkaa	75004116	-	-	-
50 ml:n Nalgene-dekantterilasi, 28 paikkaa	75004129	-	-	-
100 ml:n Nalgene-dekantterilasi, 15 paikkaa	75004118	-	-	-
250 ml:n Nalgene-dekantterilasi, 9 paikkaa	75004119	-	-	-
400 ml:n Nalgene-dekantterilasi, 8 paikkaa	75004120	-	-	-
600 ml:n Nalgene-dekantterilasi, 5 paikkaa	75004121	-	-	-
Коко 9 х 18			Telineet/paikat	
30 ml:n Nalgene-dekantterilasi, 32 paikkaa	75004110	3 / 96	4 / 128	6 / 192
50 ml:n Nalgene-dekantterilasi, 24 paikkaa	75004128	3/72	4 / 96	6/144
100 ml:n Nalgene-dekantterilasi, 15 paikkaa	75004112	3 / 45	4 / 60	6 / 90
250 ml:n Nalgene-dekantterilasi, 8 paikkaa	75004113	3/24	4 / 32	6/48
400 ml:n Nalgene-dekantterilasi, 6 paikkaa	75004114	3/18	4 / 24	6/36
600 ml:n Nalgene-dekantterilasi, 4 paikkaa	75004115	3/12	4 / 16	6/24

Taulukko 33: Saatavilla olevat Nalgene-dekantterilasin telineet Solaris 4000 -alustoille

# Solaris 2000 I / 2000 R -alustoille

	Tuotenro	Yleinen
Koko 11 x 14 / 12 x 14		Telineet/paikat
30 ml:n Nalgene-dekantterilasi, 36 paikkaa	75004116	1/36
50 ml:n Nalgene-dekantterilasi, 28 paikkaa	75004129	1 / 28
100 ml:n Nalgene-dekantterilasi, 16 paikkaa	75004118	1 / 16
250 ml:n Nalgene-dekantterilasi, 9 paikkaa	75004119	1/9
400 ml:n Nalgene-dekantterilasi, 8 paikkaa	75004120	1/8
600 ml:n Nalgene-dekantterilasi, 5 paikkaa	75004121	1/5
Koko 9 x 18		Telineet/paikat
30 ml:n Nalgene-dekantterilasi, 32 paikkaa	75004110	-
50 ml:n Nalgene-dekantterilasi, 24 paikkaa	75004128	-
100 ml:n Nalgene-dekantterilasi, 15 paikkaa	75004112	-
250 ml:n Nalgene-dekantterilasi, 8 paikkaa	75004113	-
400 ml:n Nalgene-dekantterilasi, 6 paikkaa	75004114	-
600 ml:n Nalgene-dekantterilasi, 4 paikkaa	75004115	-

Taulukko 34: Saatavilla olevat Nalgene-dekantterilasin telineet Solaris 2000 I / 2000 R -alustoille
Tarvikkeet

# Solaris 4000 I / 4000 R -alustoille

	Tuotenro	Yleinen
		18 x 18
Koko 11 x 14 / 12 x 14		Telineet/paikat
30 ml:n Nalgene-dekantterilasi, 36 paikkaa	75004116	-
50 ml:n Nalgene-dekantterilasi, 28 paikkaa	75004129	-
100 ml:n Nalgene-dekantterilasi, 16 paikkaa	75004118	-
250 ml:n Nalgene-dekantterilasi, 9 paikkaa	75004119	-
400 ml:n Nalgene-dekantterilasi, 8 paikkaa	75004120	-
600 ml:n Nalgene-dekantterilasi, 5 paikkaa	75004121	-
Koko 9 x 18		Telineet/paikat
30 ml:n Nalgene-dekantterilasi, 32 paikkaa	75004110	2 / 64
50 ml:n Nalgene-dekantterilasi, 24 paikkaa	75004128	2 / 48
100 ml:n Nalgene-dekantterilasi, 15 paikkaa	75004112	2/30
250 ml:n Nalgene-dekantterilasi, 8 paikkaa	75004113	2 / 16
400 ml:n Nalgene-dekantterilasi, 6 paikkaa	75004114	2 / 12
600 ml:n Nalgene-dekantterilasi, 4 paikkaa	75004115	2/8

Taulukko 35: Saatavilla olevat Nalgene-dekantterilasin telineet Solaris 4000 I / 4000 R -alustoille

# 1. 2. 9. Erotussuppilon pidike

### Solaris 2000 -alustoille



	Tuotenro	Yleinen		einen Dual Stack		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Erotussuppilon yksittäispidike, 250–2 000 ml	75004125	2	З	4	2	5
Pystysuoraan asennettu erotussuppilon pidike		4	4	6	-	-

Taulukko 36: Saatavilla olevat erotussuppilon pidikkeet Solaris 2000 -alustoille

# Solaris 4000 -alustoille

	Tuotenro	Yleinen		Dual Stack
		18 x 30	36 x 24	18 x 30
Erotussuppilon yksittäispidike, 250–2 000 ml	75004125	5	7	9
Pystysuoraan asennettu erotussuppilon pidike		8	12	-

Taulukko 37: Saatavilla olevat erotussuppilon pidikkeet Solaris 4000 -alustoille

# Solaris 2000 I / 2000 R -alustoille

	Tuotenro	Yleinen
		11 x 14
Erotussuppilon yksittäispidike, 250–2 000 ml	75004125	-
Pystysuoraan asennettu erotussuppilon pidike		-

Taulukko 38: Saatavilla olevat erotussuppilon pidikkeet Solaris 2000 I / 2000 R -alustoille

#### Solaris 4000 I / 4000 R -alustoille

	Tuotenro	Yleinen
		18 x 18
Erotussuppilon yksittäispidike, 250–2 000 ml	75004125	-
Pystysuoraan asennettu erotussuppilon pidike		-

Taulukko 39: Saatavilla olevat erotussuppilon pidikkeet Solaris 4000 I / 4000 R -alustoille

Tarvikkeet

# 1. 2. 10. Tarviketaso

# Solaris 2000 -alustoille



	Tuotenro	Yleinen			Dual	Stack
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
11 x 14 / 12 x 14 Tarviketaso	75004123	1	-	-	2	-
9 x 18 Tarviketaso	75004122	-	2	3	-	4

Taulukko 40: Saatavilla olevat tarviketasot Solaris 2000 -alustoille

### Solaris 4000 -alustoille

	Tuotenro	Yleinen		Dual Stack
		18 x 30	36 x 24	18 x 30
11 x 14 / 12 x 14 Tarviketaso	75004123	-	-	-
9 x 18 Tarviketaso	75004122	3	6	6

Taulukko 41: Saatavilla olevat tarviketasot Solaris 4000 -alustoille

# Solaris 2000 I / 2000 R -alustoille

	Tuotenro	Yleinen
		11 x 14
11 x 14 / 12 x 14 Tarviketaso	75004123	1
9 x 18 Tarviketaso	75004122	-

Taulukko 42: Saatavilla olevat tarviketasot Solaris 2000 I / 2000 R -alustoille

# Solaris 4000 I / 4000 R -alustoille

	Tuotenro	Yleinen
		18 x 18
11 x 14 / 12 x 14 Tarviketaso	75004123	-
9 x 18 Tarviketaso	75004122	2

Taulukko 43: Saatavilla olevat tarviketasot Solaris 4000 I / 4000 R -alustoille

#### Tarviketason lisätarvikkeet

	Tuotenro	Yleinen
O-renkaiden varaosasarja (tarviketaso)	75004132	-

Taulukko 44: Saatavilla olevat tarviketason lisätarvikkeet

# 1. 2. 11. Liimakiinnitteiset matot

Yleisiä huomautuksia liimakiinnitteisten mattojen valinnasta:

- 280 x 170 mm:n maton koko on optimoitu kokojen 11 x 14 ja 12 x 14 alustoille.
- 200 x 200 mm:n maton koko on optimoitu suuremmille kokojen 18 x 18, 18 x 24, 30 x 18 ja 36 x 24 alustoille.
- Kaikki matot voidaan leikata haluttuun kokoon.

# Solaris 2000 -alustoille

	Tuotenro	Yleinen		n Dual Stac Universa		Stack ersal
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
200 x 200 mm:n tiukasti kiinnittyvä matto, 2 kpl:n pakkaus	75004126	1	4	6	2	8
280 x 170 mm:n tiukasti kiinnittyvä matto, 2 kpl:n pakkaus	75004127	2	2	4	4	4
200 x 200 mm:n löyhästi kiinnittyvä matto, 2 kpl:n pakkaus	75004111	1	4	6	2	8
280 x 170 mm:n löyhästi kiinnittyvä matto, 2 kpl:n pakkaus	75004117	2	2	4	4	4

Matot myydään kahden kappaleen pakkauksina. Alustalle sopivien mattojen määrät eivät koske tätä pakkauskokoa vaan yhtä mattoa.

Taulukko 45: Saatavilla olevat liimakiinnitteiset matot Solaris 2000 -alustoille

# Solaris 4000 -alustoille

	Tuotenro	Yleinen		Dual Stack
		18 x 30	36 x 24	18 x 30
200 x 200 mm:n tiukasti kiinnittyvä matto, 2 kpl:n pakkaus	75004126	6	12	12
280 x 170 mm:n tiukasti kiinnittyvä matto, 2 kpl:n pakkaus	75004127	6	9	12
200 x 200 mm:n löyhästi kiinnittyvä matto, 2 kpl:n pakkaus	75004111	6	12	12
280 x 170 mm:n löyhästi kiinnittyvä matto, 2 kpl:n pakkaus	75004117	6	9	12

Matot myydään kahden kappaleen pakkauksina. Alustalle sopivien mattojen määrät eivät koske tätä pakkauskokoa vaan yhtä mattoa.

Taulukko 46: Saatavilla olevat liimakiinnitteiset matot Solaris 4000 -alustoille

# Solaris 2000 I / 2000 R -alustoille

	Tuotenro	Yleinen
		11 x 14
200 x 200 mm:n tiukasti kiinnittyvä matto, 2 kpl:n pakkaus	75004126	1
280 x 170 mm:n tiukasti kiinnittyvä matto, 2 kpl:n pakkaus	75004127	2
200 x 200 mm:n löyhästi kiinnittyvä matto, 2 kpl:n pakkaus	75004111	1
280 x 170 mm:n löyhästi kiinnittyvä matto, 2 kpl:n pakkaus	75004117	2

Matot myydään kahden kappaleen pakkauksina. Alustalle sopivien mattojen määrät eivät koske tätä pakkauskokoa vaan yhtä mattoa.

Taulukko 47: Saatavilla olevat liimakiinnitteiset matot Solaris 2000 I / 2000 R -alustoille

# Solaris 4000 I / 4000 R -alustoille

	Tuotenro	Yleinen
		18 x 18
200 x 200 mm:n tiukasti kiinnittyvä matto, 2 kpl:n pakkaus	75004126	4
280 x 170 mm:n tiukasti kiinnittyvä matto, 2 kpl:n pakkaus	75004127	2
200 x 200 mm:n löyhästi kiinnittyvä matto, 2 kpl:n pakkaus	75004111	4
280 x 170 mm:n löyhästi kiinnittyvä matto, 2 kpl:n pakkaus	75004117	2

Matot myydään kahden kappaleen pakkauksina. Alustalle sopivien mattojen määrät eivät koske tätä pakkauskokoa vaan yhtä mattoa.

Taulukko 48: Saatavilla olevat liimakiinnitteiset matot Solaris 4000 I / 4000 R -alustoille

Tarvikkeet

Con

# 1. 2. 12. Yleiset lisätarvikkeet

	Tuotenro	
Liimakiinnitteisen maton kiinnitystela	75004124	
Ruuvitaltta lisätarvikkeille	75004131	
Kaasunjakoputkisto (Solaris 2000 I / 2000 R)	SK2000-8GM	
Kaasunjakoputkisto (Solaris 4000 I / 4000 R)	SK4000-8GM	
Puristimen varaosasarja (Ruuveja)	SK0010	
Alustan varaosasarja alustoille Solaris 2000, 2000 I/R ja 4000 I/R (alustan ruuvit, työkalu)	SK0100	
Alustan varaosasarja SK4000:lle (Alustan ruuvit, työkalu)	SK0101	

Taulukko 49: Saatavilla olevat yleiset lisätarvikkeet

# 1. 3. Direktiivit ja standardit

Alue	Direktiivi	Standardit		
Eurooppa	2006/42/EY	EN 61010-1		
	Konedirektiivi	IEC 61010-2-051		
	2011/65/EU RoHS	EN 61326-1 Luokka B		
	Direktiivi tiettyjen vaarallisten	EN ISO 14971		
	aineiden käytönrajoittamisesta sähkö- jaelektroniikkalaitteissa	EN ISO 9001		
	Suojelukohteet:	Ainoastaan alustoille: Solaris		
	2014/35/EU Alhainen jännite	2000 17 2000 R7 4000 17 4000 B' IEC 61010-2-010		
	2014/30/EY Sähkömagneettinen yhteensopivuus (EMC)			
Pohjois Amerikka		ANSI/UL 61010-1		
		IEC 61010-2-051		
		IEC 61326-1 Luokka B		
		CFR 47 FCC 15 EMC		
		EN ISO 14971		
		EN ISO 9001		
		Ainoastaan alustoille: Solaris 2000 I / 2000 R / 4000 I / 4000 R: IEC 61010-2-010		
Japani		IEC 61010-1		
		IEC 61010-2-051		
		IEC 61326-1 Luokka B		
		EN ISO 14971		
		EN ISO 9001		
		Ainoastaan alustoille: Solaris 2000 I / 2000 R / 4000 I /		
		4000 R: IEC 61010-2-010		

Taulukko 50: Direktiivit ja standardit

HUOMIO: Tämä laite on testattu ja sen on havaittu noudattavan rajoja luokan A digitaalisille laitteille, FCC-sääntöjen kohdan 15 mukaan. Nämä rajat on kehitetty kohtuulliseen suojaamiseen haitallista interferenssiä vastaan kun laitetta käytetään kaupallisessa ympäristössä. Tämä laite tuottaa, käyttää ja voi aiheuttaa radiofrekvenssienergiaa ja voi, jos se ei ole asianmukaisesti asennettu ja käytetty käyttöoppaan mukaan, aiheuttaa haitallista interferenssiä radioyhteyksille. Tämän laitteen käyttö asuinalueella aiheuttaa mahdollista haitallista interferenssiä, jossa tapauksessa käyttäjä on velvollinen korjaamaan interferenssin omalla kustannuksellaan.

# 1.4. Tuotekatsaus

# 1.4.1. Solaris 2000



① Alustan kiinnityslevy; ② Graafinen käyttöliittymä; ③ Virtakatkaisin; ④ Verkkovirtaliitäntä;
 ⑤ Sulake; ⑥ Ethernet-portti; ⑦ USB-portti

Kuva 1: Yleiskuva Solaris 2000

## 1.4.2. Solaris 4000



① Alustan kiinnityslevy; ② Graafinen käyttöliittymä; ③ Virtakatkaisin; ④ Verkkovirtaliitäntä;
 ⑤ Sulake (uudelleenasetettava); ⑥ Ethernet-portti; ⑦ USB-portti

Kuva 2: Yleiskuva Solaris 4000

Tuotekatsaus

# 1. 4. 3. Solaris 2000 I / 2000 R



① Alustan kiinnityslevy; ② Graafinen käyttöliittymä; ③ Virtakatkaisin; ④ Käyttöportti;
⑤ Verkkovirtaliitäntä; ⑥ Sulake (uudelleenasetettava); ⑦ Ethernet-portti; ⑧ USB-portti;
⑨ Kansi

Kuva 3: Solaris 2000 I / 2000 R - Yleiskatsaus

# 1.4.4. Solaris 4000 I / 4000 R



① Alustan kiinnityslevy; ② Graafinen käyttöliittymä; ③ Virtakatkaisin; ④ Käyttöportti;
⑤ Verkkovirtaliitäntä; ⑥ Sulake (uudelleenasetettava); ⑦ Ethernet-portti; ⑧ USB-portti;
⑨ Kansi

Kuva 4: Solaris 4000 I / 4000 R - Yleiskatsaus

# 1.4.5. Liitännät

### Avoimet ravistimet



1) Verkkovirtaliitäntä; 2) Sulake; 3) Ethernet-portti

Kuva 5: Takanäkymä (vas.: Solaris 4000, oik.: Solaris 2000)

# Inkuboidut ja jäähdytetyt ravistimet



① Verkkovirtaliitäntä; ② Sulake; ③ Ethernet-portti; ④ Tyhjennysputki
 Kuva 6: Takanäkymä (vas.: Solaris 2000 I / 2000 R, oik.: Solaris 4000 I / 4000 R)

# <u>Verkkovirtaliitäntä</u>

Ravistin tarvitsee teknisten tietojen mukaisen teholähteen. Virtakaapelit tulevat mukana toimituksessa.

Tuotenro	Ravistin	Tekniset tiedot
SK2000	Solaris 2000	100–240 V ±10 %, 50 / 60 Hz
SK4000	Solaris 4000	100–240 V ±10 %, 50 / 60 Hz
SK2001	Solaris 2000 I	100–120, 200–240 V ±10 %, 50 / 60 Hz
SK4001	Solaris 4000 I	100–120, 200–240 V ±10 %, 50 / 60 Hz
SK2002	Solaris 2000 R	100–240 V ±10 %, 50 / 60 Hz
SK4002	Solaris 4000 R	100–240 V ±10 %, 50 / 60 Hz

Kuva 7: Ravistimien teholähteiden tekniset tiedot

VAROITUS Väärä verkkoliitäntäpistoke tai pistotulppa aiheuttaa vaurion. Varmista, että ravistin liitetään ainoastaan sellaisiin pistorasioihin, jotka on maadoitettu asianmukaisesti voimassa olevien sähköturvallisuusstandardien mukaisesti. Älä käytä ravistinta viallisella tai liian alhaisen luokituksen omaavalla virtajohdolla.

Pääkatkaisijan on oltava vapaasti saavutettavissa kaikkina aikoina.

Estääksesi sähköiskun vaaran huolehdi, että ympäröivät pinnat ovat kuivat. Jos nesteitä pääsee vahingossa läikkymään tai roiskumaan, irrota ravistin teholähteestä, puhdista läikkyneet nesteet ja torju kaikki terveys- tai biologiset vaarat ennen kuin jatkat työskentelyä.

Kytke ravistin irti virtalähteestä kun se ei ole käytössä.

#### <u>Sulake</u>

Sulake voidaan kytkeä takaisin päälle, jos se on lauennut ylivirran takia. Ravistin ei toimi, ennen kuin sulake on työnnetty takaisin oikeaan asentoon.

#### Ethernet

Ravistimessa on RJ45 Ethernet -portti, jota voidaan käyttää lähiverkkoon (LAN) liittymiseen. Käytä RJ45 Ethernet -portissa ainoastaan laitteita, jotka täyttävät standardin IEC 60950-1 vaatimukset. RJ45 Ethernet -portti on valmis käyttöön jatkossa, kun sopiva ohjelmistopäivitys on saatavilla.

#### <u>Yhteysportit</u>

Lämpötilaohjatuissa ravistinmalleissa on kaksi sivuaukkoa, joita kutsutaan yhteysporteiksi. Nämä portit on yleensä tulpattu korkeilla. Voit irrottaa korkin ja ottaa käyttöön muita laitteita, kuten ylimääräisen lämpötila-anturin, tai syöttää kaapelin tai letkun, esimerkiksi kaasunjakoputkiston putket.

# <u>USB</u>

Ravistimessa on kaksi USB-A 2.0 -porttia, joihin sopivat kaupallisesti saatavilla olevat USB-muistit. Käytä vain laitteita jotka täyttävät standardissa IEC 60950-1 USB-porteille asetetut vaatimukset.

#### Sisäinen ylikuumenemissuojaus

Lämpötilaohjatuissa ravistimissa on lämmityselementtien viereen asennettu manuaalisesti nollattava termostaatti. Tämä lämpölaite sammuttaa lämmittimet, jos ilmaa kierrättävät puhaltimet vikaantuvat, jotta lämpötila ei nouse liian korkeaksi ravistimen kotelon sisällä.

Manuaalinen termostaatti tulee nollata käsin, jotta ravistin saadaan taastoimintakuntoon. Tämän voi tehdä ainoastaan Thermo Fisher Scientific -asiakaspalvelu.

# 2. Kuljetus ja asennus



On sinun vastuullasi varmistaa että ravistin on asennettu asianmukaisesti.

#### HUOMIO

Toimituspakkaus pitää tutkia sen saapuessa. Tutki tarkoin mahdolliset toimitusvauriot ennen pakkauksen purkamista. Jos vaurioita havaitaan, toimittajan tulee eritellä ne ja allekirjoittaa vaurioiden löytyminen teidän kappaleeseenne toimituskuitista.

Avaa pakkaus varovasti ja tarkista, että kaikki osat ("Taulukko 51: Tuotteet") ovat mukana toimituksessa ennen kuin hävität pakkausmateriaalit. Purkamisen jälkeen, jos vaurioita löytyy, ilmoita niistä kuljetusyhtiölle ja vaadi vauriotarkastusta.

Tärkeää: Jos et pyydä vauriotarkastusta muutaman päivän sisällä toimituksesta, kuljetusyhtiö vapautuu vaurion vastuusta. Sinun on pyydettävä vauriotarkastus.

# 2.1. Purku

Käytä pakkauslistaa, kun purat, todentaaksesi että koko laite on vastaanotettu. Älä heitä pakkausmateriaaleja pois ennen kuin kaiken on todettu olevan mukana.

# Tuotteet

Esine	Määrä
Ravistin	1
Virtajohto	1
Yleisalusta	1*
Alustan ruuvit	
- Solaris 4000	4*
- muut mallit	3*
Painettu englanninkielinen	1
käyttöopas	
Käyttöohjeet USB:llä	1
Lukitustyökalu (T-kahvainen avain)	1
* 2000 I-, 2000 R-, 4000 I- ja 4000 R -ravis	timissa on ravistimeen
valmiiksi asennettu alusta. Alustaa tai ruuve	ja ei ole saatavana erillisinä
osina.	

Taulukko 51: Tuotteet

Jos jotakin puuttuu, ota yhteys Thermo Fisher Scientificiin.

# 2.2. Sijoituspaikka

NOUDATA VAROVAISUUTTA Suojaus voi vaarantua alentuneen stabiliteetin myötä muovin altistuttua ultraviolettisäteilylle. Älä altista ravistinta ja muoviapuvälineitä suoralle auringonvalolle tai muille ultraviolettisäteilyn lähteille.

Aseta ravistin tasaiselle pöydälle tai penkille, joka pystyy kantamaan ravistimen ja mahdolliset apuvälineet ja näytteet sen ollessa käytössä. Aseta ravistin lähelle virtalähdettä joka vastaa arvolevyn vaatimuksia. Jätä tyhjää tilaa laitteen ympärille ilmanvaihdolle, apuvälineiden kiinnitystä varten ja käyttäjälle.

Ota huomioon seuraavat vaatimukset asentaessasi laitetta:

- Avoin ravistin tarvitsee kaikille sivuilleen vapaata tilaa 8 cm (3 tuumaa), jotta liikkuville osille jää riittävä välys ja jotta vältytään puristumisvahingoilta ja viereisten laitteiden vaurioitumiselta. Avoimen ravistimen alusta voi mennä päällekkäin kotelon kanssa.
- Ravistin aiheuttaa tärinää. Älä säilytä herkkiä laitteita tai vaarallisia esineitä tai aineita turvavyöhykkeellä.



Kuva 8: Asennus 8 cm:n (3 tuuman) etäisyydelle avoimesta ravistimesta

Lämpötilaohjattujen ravistimien osalta tulee noudattaa seuraavia tilavaatimuksia, jotta tuuletus toimisi oikein ja suorituskyky olisi paras mahdollinen:

- Inkuboidut Solaris 2000 I ja 4000 I ravistimet vaativat 8 cm:n (3 tuuman) vapaan tilan kotelon kaikilla neljällä sivulla Kuva 9 vasemmalla puolella näkyvällä tavalla.
- Jäähdytettyjä Solaris 2000 R ja 4000 R -ravistimia voidaan käyttää niin, että kotelon kaikilla neljällä sivulla on 8 cm:n (3 tuuman) vapaa tila, jos niiden kannen yläpuolella on vapaata ilmatilaa Kuva 9 vasemmalla puolella näkyvällä tavalla.

#### Sijoituspaikka

 Jos jäähdytetyn Solaris 2000 R - ja 4000 R -ravistimen yläpuolella oleva ilmatila on tukkeutunut, kotelon ympärillä oleva vapaa tila tulee kasvattaa 30 cm:iin (12 tuumaan) kaikilla neljällä sivulla Kuva 9 oikealla puolella esitetyllä tavalla.



**Kuva 9:** Määritetyt vapaat tilat inkuboiduille (vasemmalla) ja jäähdytetyille (oikealla) ravistimille. Jäähdytettyjen vapaa ilmatila vasemman kuvan mukaisesti.

VAROITUS Vaaralliset aineet on pidettävä pois tältä alueelta sekoitettaessa.

Kantavan pinnan täytyy:

- » olla puhdas,
- » olla vakaa, tukeva, jäykkä ja tärinätön,
- » soveltua ravistimen asentamiseen vaakasuunnassa,
- » kantaa ravistimn paino.
- Ravistinta tulee käyttää vain sisätiloissa.
- Ravistinta ei tule altistaa kuumuudelle ja voimakkaalle auringonvalolle.
- Laitteen sijainnin tulee olla aina hyvin tuuletettu.
- Pääkatkaisijan on oltava vapaasti saavutettavissa kaikkina aikoina.

# 2.3. Kuljetus

NOUDATA VAROVAISUUTTA Fyysinen vahinko, joka johtuu laitteen pudottamisesta. Nosta aina ravistinta molemmilta puolilta. Älä koskaan nosta ravistinta etupaneelista tai asennetusta alustasta. Älä nosta lämpötilaohjattua ravistinta läpinäkyvän kannen kahvasta.

NOUDATA VAROVAISUUTTA Kotelon sisälle tiivistyvä vesi vaikuttaa haitallisesti ravistimen turvalliseen käyttöön. Näin tapahtuu, kun ravistin siirretään viileämmästä lämpimämpään paikkaan, jos ilma on kosteampaa, esimerkiksi varastoalueelta tai kuljetusautosta laboratoriotilaan. Anna laitteen kuivua vähintään 2 h ennen käyttöä.

HUOMIO Poista aina avoimen ravistimen alusta, sen kuormitus ja kaikki lisätarvikkeet ennen ravistimen siirtämistä. Jos et poista alustaa, saatat vahingoittaa alustan kiinnityslevyä tai sekoitusmekanismia. Lämpötilaohjatuissa ravistimissa riittää, että kuormitus poistetaan alustalta.

**HUOMIO** Yksikkö voi vaurioitua iskusta. Kuljeta laitetta pystyasennossa ja jos mahdollista, alkuperäispakkauksessa.

# Ravistimen käsittely

Huolehdin ravistimen käsittelyn yhteydessä siitä, että:

- nostat ravistinta kummaltakin puolelta mutta et koskaan edestä etkä takaa
- annat ravistimen niin monen henkilön kannettavaksi kuin ravistimen paino edellyttää (katso kohtaa "Tekniset tiedot" sivulla 11) mutta aina kuitenkin vähintään kahdelle henkilölle





Kuva 10: Miten ravistinta kannetaan

VAROITUS Nosta aina ravistinta molemmilta puolilta. Älä nosta ravistinta edestä tai takaa. Ravistimet ovat raskaita (katso kohtaa "Tekniset tiedot" sivulla 11). Anna ravistin vähintään kahden henkilön nostettavaksi ja kannettavaksi.

**NOUDATA VAROVAISUUTTA** Älä nosta lämpötilaohjattua ravistinta läpinäkyvän kannen kahvasta. Kahva voi murtua ja ravistin voi pudota maahan ja aiheuttaa vakavan henkilövamman.

# 2.4. Vaaitus

Ravistinta ei voi vaaittaa yksinään. Ravistin vaatii vaaitetun tukipinnan kunnolliseen asennukseen.

HUOMIO Älä pane levyjä tai muita litteitä esineitä ravistimen jalkojen ja niiden alustan väliin ravistimen vaakasuoraan saamiseksi.

# 2.5. Verkkovirtaliitäntä

Ravistin tarvitsee teknisten tietojen mukaisen teholähteen. Virtakaapelit tulevat mukana toimituksessa. Katso lisätietoja Kuva 7 sivulla 47.

VAROITUS Vahinko väärän verkkoliitäntäpistokkeen tai virransaantipistokkeen käytöstä. Pidä huoli että ravistin on kytketty vain kunnolla maadoitettuihin pistorasioihin. Älä käytä ravistinta viallisella tai liian alhaisen luokituksen omaavalla virtajohdolla.

HUOMIO Sähkömagneettinen säteily voi aiheuttaa häiriöitä näytössä. Tämä ei vahingoita laitetta, rajoita sen toimintaa tai muuta laitetta. Älä tuo mobiililaitteita, kuten matkapuhelimia, laitteen lähelle, jotta ne eivät aiheuta sähkömagneettisesta säteilystä aiheutuvaa häiriötä. Älä käytä laitetta samassa virtapiirissä muiden paljon virtaa kuluttavien laitteiden kanssa. Älä käytä useita laitteita samassa jatkojohdossa.

Yhdistääksesi ravistimen virrantuloon tee seuraavasti:

- 1. Käännä oikealla sivulla sijaitseva virtakytkin pois päältä.
- 2. Varmista, että virtajohto vastaa oman maasi standardeja.
- 3. Varmista, että jännite ja taajuus ovat samat kuin laitekilvessä.

Pääkatkaisijan on oltava vapaasti saavutettavissa kaikkina aikoina.

Estääksesi sähköiskun vaaran huolehdi, että ympäröivät pinnat ovat kuivat. Jos esiintyy tahatonta nesteiden läikkymistä tai roiskumista, kytke ravistin irti virtalähteestä, puhdista läikkyneet nesteet ennen kuin jatkat työskentelyä.

Kytke ravistin irti virtalähteestä kun se ei ole käytössä.

# 2.6. Ensimmäinen käynnistys

Käy seuraavat vaiheet läpi ennen kuin käytät laitetta:

1. Pane virta päälle, thermo scientific- logo näyttäytyy. Napauta Start Setup.



Kuva 11: Alkukäynnistyksen kehote

2. Valitse kieli Kielinäytössä. Napauta Seuraava.



Kuva 12: Ensimmäinen käynnistys - kieli

 Voit halutessasi antaa yksikölle nimen valintaikkunassa Yksikön nimi. Napauta Seuraava.

	Unit N	ame	
Unit Name:	Tap to enter		
E	Back	Next	

Kuva 13: Ensimmäinen käynnistys - laitteen nimi

 Voit halutessasi syöttää kaupungin ja maan tekstikenttään. Kun olet kirjoittanut kolme merkkiä, Kaupunki ja maa -tekstikentässä näkyy vastaavien ehdotusten luettelo. Voit halutessasi valita jonkin ehdotuksen ja napauttaa sen jälkeen Seuraava.

Set Re	gion						
Enter the city and co	untry closest to your time zone						
City, Country: Tap to enter							
ack	Next						
	Set Re Enter the city and co Tap to ente	Set Region Enter the city and country closest to your time zone Tap to enter					

Kuva 14: Ensimmäinen käynnistys - alue

5. Valitse haluttu päiväysformaatti. Napauta Seuraava.

Set Date Dis	olay Format							
MMN	1/DD/YYYY							
O DD/N	O DD/MMM/YYYY							
O YYYY	/MMM/DD							
Back	Next							

Kuva 15: Ensimmäinen käynnistys – aseta päiväysformaatti

6. Valitse nykyinen päivämäärä kääntämällä valintakiekkoa. Napauta Seuraava.



Kuva 16: Ensimmäinen käynnistys - aseta päiväys

7. Valitse haluamasi ajan muoto ja määritä nykyinen aika kääntämällä valintakiekkoa. Napauta **Seuraava**.



Kuva 17: Ensimmäinen käynnistys – aseta aika

8. Vain inkuboidut ja jäähdytetyt ravistimet: Valitse haluamasi lämpötilan yksikkö. Napauta Seuraava.

Temperature Units								
۲	°C							
0	°F							
Back	Next							

Kuva 18: Alkukäynnistys – valitse lämpötilan yksikkö

 Vain inkuboidut ja jäähdytetyt ravistimet: Aseta korkean lämpötilan hälytyksen ja hätäpysäytyksen kynnysarvot napsauttamalla sinistä nuolta. Napauta Seuraava.



Kuva 19: Alkukäynnistys – aseta korkean ja matalan lämpötilan hälytyskynnykset

10. Toista vaiheet matalan lämpötilan hälytyskynnysten osalta. Jatka napauttamalla **Seuraava**.

11. Jos haluat jatkaa asettamatta järjestelmänvalvojan salasanaa, jätä Avoin tila valituksi ja jatka valitsemalla Seuraava.

Jos haluat määrittää järjestelmänvalvojan salasanan nyt, napauta Suojattu tila.

	Acces	s Mode
Op Sele	en Mode ect to allow full access t	o all functionality for all users.
O Sel Sel cha (All	cure Mode ect to require users to en nges to stored paramet users still able to use pro	nter the admin passcode to make ers, programs, and settings. grams and set new run parameters,
	Back	Next

Kuva 20: Alkukäynnistys – Valitse käyttötila

12. Kirjoita näyttöön tulevaan salasanakehotteeseenjärjestelmänvalvojan salasana näppäimistöltä ja napauta **Seuraava**.



Kuva 21: Alkukäynnistys – Valitse käyttötila

- 13. Kirjoita järjestelmänvalvojan salasana seuraavaan näyttöön tulevaan salasanakehotteeseen.
- 14. Anna järjestelmänvalvojan salasana uudelleen vielä kerran, kun näyttöön tulee kolmas salasanakehote.

	Admin Passcode					Admin Passcode								
If you	do not wis fields	h to change below blanl	the pa and ta	sscode and the Ne	at this ti ext butto	me, leav on.	e the		F	Please re-en	ter the n	ew passco	de.	
	*	*	*		*	*			*	*	*	*	*	
		7	8	9						7	8	9		
		4	5	6						4	5	6		
		1	2	3						1	2	3		
		0	Cle	ear						0	Cle	ar		

Kuva 22: Alkukäynnistys – Anna uusi järjestelmänvalvojan salasana ja vahvista se

15. Jatka valitsemalla näkyviin tulevassa vahvistusnäytössä Seuraava.

Access Mode			
New passcode create successfully. Please keep the new passcode in a convenient and safe location as you will need it to change settings.			
If the passcode is forgotten you will need to contact Customer Service to reset.			
Back	Next		

Kuva 23: Alkukäynnistys – Järjestelmänvalvojan salasanan muutoksen vahvistus

 Näyttöön tulee ikkuna, jossa on ravistelijan yleiset fyysiset asennusohjeet. Napauta Seuraava.

Installation Instructions			
Shakers must be installed on a stable and level surface with sufficient air circulation. Please reference the product manual for instruction on the proper installation of this unit.			
Back	Next		

Kuva 24: Ensimmäinen käynnistys - asennusohje

17. Näyttöön tulee Asennus valmis -ikkuna. Lopeta napauttamalla Seuraava.

Setup Complete				
Setup is complete.				
We recommend calibrating the unit at the ideal temperature for your application. Would you like to calibrate the temperature now? Calibration can always be accessed under the main settings screen.				
Calibrate	Finish			

Kuva 25: Alkukäynnistys – kalibroi tai viimeistele asetukset

HUOMIO Inkuboiduissa ja jäähdytetyissä ravistimissa voit napauttaa **Kalibroi** ja suorittaa toimenpiteet, jotka on lueteltu luvussa "4. 6. Lämpötilan kalibrointi" sivulla 126.

# 2.7. Varastointi



ja jos tarpeen, desinfioi tai dekontaminoi koko systeemi. Älä jätä ravistinta ja apuvälineitä määrittelemättömään NOUDATA kontaminaatiotilaan. Jos et ole varma prosessista, ota yhteys VAROVAI-Thermo Fisher Scientificin asiakaspalveluun ("Puhdistus" sivulla SUUTTA 123, "Desinfiointi" sivulla 124 ja "Dekontaminaatio" sivulla 124).

Kun otat ravistimen ja apuvälineet pois käytöstä, puhdista

- Ennen kuin varastoit ravistimen ja apuvälineet, ne on puhdistettava, ja jos on tarpeen, desinfioitava ja dekontaminoitava.
- Ravistimen ja apuvälineiden on oltava täysin kuivat ennen varastointia.
- Pidä ravistin puhtaassa pölvttömässä paikassa.
- Pidä ravistin jaloillaan.
- Älä varastoi ravistinta suorassa auringonvalossa.

# 2.8. Toimitus



NOUDATA

VAROVAI-

SUUTTA

Ennen kuljetusta ravistin ja apuvälineet on puhdistettava ja jos tarpeen, desinfioitava ja dekontaminoitava koko järjestelmä. Älä jätä ravistinta ja apuvälineitä määrittelemättömään kontaminaatiotilaan. Jos et ole varma prosessista, ota yhteys Thermo Fisher Scientificin asiakaspalveluun ("Puhdistus" sivulla 123, "Desinfiointi" sivulla 124 ja "Dekontaminaatio" sivulla 124).

Ennen ravistimen kuljetusta:

- Ravistimen tulee olla puhdas ja dekontaminoitu.
- Dekontaminaatio on varmennettava dekontaminaatiotodistuksella. Dekontaminaatiotodistuksia saa Thermo Fisher Scientificin asiakaspalvelusta.

# 3. Käyttö

# 3.1. Virta päälle / pois

Paina virtakytkin ravistimen oikealla puolella päälle (I) tai pois (0).

Kosketusnäytöllä on Thermo Scientific -logo kun se menee päälle.

Kun se on valmis, kosketusnäyttö näyttää ravistimen tämänhetkisen tilan.

# 3. 2. Graafinen käyttöliittymä

Aloitusnäyttö on ravistimen graafisen käyttöliittymän (GUI) oletusnäyttö. Ravistimen käyttö aloitetaan siitä.

Aloitusnäytössä voit:

- määrittää käytön perusparametrit, kuten ravistusnopeuden, ajan ja lämpötilan (vain lämpötilaohjatut ravistimet)
- käynnistää ja pysäyttää ravistimen
- tarkastella tilatietoja ja käsitellä hälytyksiä ja hälytysilmoituksia
- siirtyä muihin näyttöihin, joissa on tilatietoja ja asetusvaihtoehtoja

Näytön sisältö on hieman erilainen avoimissa ja lämpötilaohjatuissa versioissa.

Avoimen ravistimen aloitusnäyttö on Kuva 26 esimerkin kaltainen.



Kuva 26: Graafisen käyttöliittymän aloitusnäyttö avoimessa ravistimessa

#### Graafinen käyttöliittymä

Lämpötilaohjatun ravistimen aloitusnäytössä on lämpötilan lisäsäätimet, jotka näkyvät Kuva 27.



Kuva 27: Graafisen käyttöliittymän aloitusnäyttö lämpötilaohjatussa ravistimessa

# 3. 2. 1. Ravistimen käytön perusparametrien määritys

Aloitusnäytöstä pääset ravistimen tärkeimmät käyttöparametreihin yhdellä napautuksella.

#### Määritä nopeus



Kuva 28: Nopeuden asetuspiste

- Määritä haluttu nopeus napauttamalla valintakiekon kunkin valitsimen ylä- tai alapuolella olevaa nuolta. Vaihtoehtoisesti voit muuttaa haluttua nopeutta kääntämällä kutakin valintakielen valitsinta erikseen.
- 3. Napauta **Tallenna**.

HUOMIO Jos näyttöön tulee ponnahdusikkuna, jossa ilmoitetaan, että nopeuden asetusarvo on alueen ulkopuolella, asetuksesi on ravistimen tuetun toimintanopeusalueen ulkopuolella, ks. kohta "Tekniset tiedot" sivulla 11. Korjaa asetus ja jatka.

 Kuittaa Asetuksesi on tallennettu -vahvistusvalintaikkuna, joka tulee näkyviin päänäyttöön palaamista varten.

#### Ratalaskuri: Laske arvioitu nopeus vanhojen laiteasetusten perusteella

Nopeuden asetusarvo -valintaikkunasta voit avata **Orbit Calculator** -ratalaskimen napauttamalla kohtaa **Laske nopeus:** Ratalaskimesta on hyötyä, kun siirryt erilaisen mekaanisen rakenteen omaavasta ravistimesta johonkin tässä käyttöoppaassa kuvattuun Solaris-sarjan ravistimeen. Tämä laskin antaa radan kokoeron perusteella karkean arvion samanlaisten tulosten saamiseksi tarvittavasta asetusnopeudesta.

HUOMIO Älä käytä Orbit Calculator -ratalaskimella saatuja tuloksia sellaisenaan tuotantonäytteissä, vaan vahvista tulokset suorittamalla sarja testejä. Solut voivat kasvaa nopeammin tai hitaammin, ilmaista eri proteiineja tai olla vaurioituneita (tai kuolla) leikkausjännitysten mukaan, kun fysiikka muuttuu siirtyessäsi erilaiselle radalle.

1. Valitse mittayksikkö millimetreissä tai tuumissa.



Kuva 29: Ratalaskuri

 Valitse edeltävän radan koko. Napauta Laske. Seuraava ikkuna ilmestyy näkyviin nopeutta laskettaessa.



Kuva 30: Rata laskettu

- 3. Napauta **OK**.
- 4. Voit laskea viitteellisen nopeuden räätälöidylle radalle. Jos haluat laskea mukautetun radan, valitse Muu.
- 5. Syötä radan koko.



Kuva 31: Laske räätälöidyn koon rata

6. Napauta Laske.

Laske nopeus-ikkuna ilmestyy näyttäen nopeuden annetulle koolle.

# <u>Määritä käyntiaika</u>

Voit käyttää ravistinta jatkuvassa tai ajastetussa tilassa. Jatkuvassa tilassa ravistin pysäytetään manuaalisesti käyttäjän harkinnan mukaan. Ajastetussa tilassa ravistimen käyttölaite pysähtyy automaattisesti, kun ajastimeen määritetty aika on kulunut loppuun. Saat ajastimen näkyviin seuraavasti:

- kulunut aika: kuinka kauan ravistin on ollut käynnissä Käynnistys-painikkeen painamisesta lähtien tai
- jäljellä oleva aika: kuinka kauan ravistin pysyy käynnissä, kun ajastimen aika on kulunut loppuun

- 1. Napauttamalla **Aika** -kenttää (⑤ Kuva 26 ja Kuva 27) voit avata Aikatila-näytön, joka näkyy Kuva 32.
- 2. Valitse **Ajastettu**.



Kuva 32: Valitse Aikaatila

- 3. Syötä tunnit ja minuutit Aseta aika -laatikoihin.
- 4. Napauta Tallenna.
- 5. Voit palata päänäyttöön kuittaamalla näyttöön tulevan **Asetuksesi on tallennettu** -vahvistusvalintaikkunan.

# Lämpötilan määritys

Lämpötilaohjatuissa malleissa **Lämpötilan asetusarvo** -kentän avulla voit määrittää käyttötarkoitukseesi sopivan lämpötilan milloin tahansa.

- Määritä esiasetettu oletuslämpötila, jonka haluat ravistimen ylläpitävän, napauttamalla valintakiekon kunkin valitsimen ylä- tai alapuolella olevia nuolia. Vaihtoehtoisesti voit muuttaa haluttua lämpötilaa kääntämällä kutakin valintakielen valitsinta erikseen.



Kuva 33: Määritä lämpötilan asetusarvo: kalibrointilukeman sisältävä näyttö

3. Napauta Tallenna.

HUOMIO Jos näyttöön tulee ponnahdusikkuna, jossa ilmoitetaan, että lämpötilan asetusarvo on alueen ulkopuolella, asetuksesi on ravistimen tuetun toimintalämpötilaalueen ulkopuolella, ks. kohta "Tekniset tiedot" sivulla 65. Korjaa asetus ja jatka.

HUOMIO Jos näyttöön tulee ponnahdusikkuna, jossa ilmoitetaan, että valittu lämpötila on ehkä alueen ulkopuolella ympäristön lämpötilan johdosta, ympäristön lämpötila on ravistimen tuetun toimintalämpötila-alueen ulkopuolella, ks. kohta "Tekniset tiedot" sivulla 11. Vahvista ilmoitus ja jatka valitsemalla **OK** tai valitse toinen lämpötila.

HUOMIO Lämpötilan asetusarvo -näytössä näkyy kalibroinnin aikana säädetty kalibroitu lämpötila ja poikkeama Kuva 33 vasemmassa alakulmassa kuvatulla tavalla. Kalibrointiprosessi on kuvattu kohdassa "Lämpötilan kalibrointi" sivulla 126.

4. Voit palata päänäyttöön kuittaamalla näyttöön tulevan **Asetuksesi on** tallennettu -vahvistusvalintaikkunan.

HUOMIO Lämpötilaohjattu ravistin toimii erityisellä tavalla, kun saavutetaan asetettu lämpötila, ns. lämpötilan ohitus. Lämpötilan ohitus tarkoittaa, että määritetty kammion lämpötila ohitetaan siirtymällä ensin hieman määritettyä korkeampaan (tai matalampaan) lämpötilaan, jonka jälkeen lähestytään määritettyä lämpötilaa. Määritetyn lämpötilan lähestymisen aikana kosketusnäyttö ei näytä tätä ohitusmenettelyä. Sen sijaan kosketusnäytössä näkyy kammion lämpötilan nousu (tai lasku), kunnes määritetty lämpötila on saavutettu.

#### Ravistimen käynnistys ja pysäytys

1. Käynnistä ravistin napauttamalla Käynnistä -painiketta.

Käynnistys-painike muuttuu Pysäytys -painikkeeksi.



Kuva 34: Käynnistys- ja Pysäytys-painike

2. Pysäytä ravistin napauttamalla painiketta Pysäytys.

# 3. 2. 2. Tila

Kun ravistin on kunnossa, kosketusnäytössä näkyy vihreän sydänkuvakkeen Kunto



ja tiedot -alueella (7) Kuva 26 ja (8) Kuva 27). Napauttamalla vihreää sydänikonia saat näkyviin tilanäytön. Tilanäytössä on tietoja ravistimen toiminnasta, mukaan lukien ravisteluun käytetyt käyttötunnit ja käyttötuntien

kokonaismäärä sekä viimeisen käytön aikainen käyttötuntimäärä. Lämpötilaohjatuissa ravistimissa näkyy myös jäähdytykseen tai lämmitykseen käytettyjen tuntien lukumäärä.

Status: <b>Good</b>	•	2	×	Status: <b>Good</b>		2	×
Total Run Time		Last Run Time Ses	ssion	Total Run Time		Last Run Time Ses	sion
Hours Shaking	232hrs	Hours Shaking	60hrs	Hours Shaking	232hrs	Hours Shaking	60hrs
Hours Powered	332bre	Hours Powered	61bre	Hours Powered	332hrs	Hours Powered	61hrs
10013 FOWERED 3321113	552113	Hous roweled only	Hrs tempering	168hrs	Hrs tempering	61hrs	
			()				(i)

Kuva 35: Ravistimen tiedot: Avoin ravistin (vasen) ja Lämpötilaohjattu ravistin (oikea)

#### Varoitus

Kun hälytys laukeaa, kosketusnäytössä näkyy keltainen vierivä tekstinauha senhetkisen näytön yläreunassa. Lisäksi kuuluu äänihälytys. Keltainen tekstinauhatyyppinen hälytyspalkki poistuu näkyvistä, kun viesti on vierinyt näytöllä kahdesti. Vain keltainen kolmio ilmaisee, että ravistimessa on yksi



tai useampi hälytys. Kolmiokuvakkeessa on valkoisella reunuksella ympäröity sininen ympyrä, joka näyttää aktiivisten hälytysten lukumäärän. Kolmiokuvakkeen painaminen Tiedot ja kunto -alueella (7) Kuva 26 ja ® Kuva 27) avaa luettelon kulloinkin voimassa olevista hälytyksistä. Viimeisin hälytys näkyy laajennettuna, mikä mahdollistaa tietojen katselun, ks. Kuva 36. Voit vierittää listaa ja napauttaa haluamaasi kohdetta laajentaaksesi sen ja lukeaksesi enemmän.



Kuva 36: Hälytysluettelo

Voit valita aktiivisen hälytyksen napauttamalla hälytyskohteen vieressä olevaa valintaruutua. Voit poistaa hälytyksen luettelosta napauttamalla painiketta **Kuittaa**. Kun kaikki hälytykset on poistettu, Tila-kuvakkeeksi palautuu oletusarvoinen vihreä sydän.

# <u>Hälytys</u>

Kun hälytys annetaan, ravistin pysähtyy välittömästi estääkseen vahingot näytteille ja/



tai laitteelle itselleen. Hälytys tulee kuitata kosketusnäytössä ennen kuin voit jatkaa työskentelyä.

Kun hälytys annetaan, kosketusnäytössä näkyy punainen palkki nykyisen näytön yläosassa. Tiedot ja kunto -alueella (⑦ Kuva 26 ja ⑧ Kuva 27) näkyy ääniaaltojen ympäröimä punainen hälytyskello. Lisäksi kuultava hälytysääni soi jatkuvasti.

Punaisen hälytyspalkin alla olevassa tekstinauhassa näkyy nykyisen hälytyksen vierivä tiivistelmä. Torkku-painike tulee näkyviin näytön oikealle puolelle ja antaa mahdollisuuden hälytyksen tilapäiseen mykistämiseen. Kun hälytystä ei selvitetä torkkujakson aikana, kuultava hälytys palaa. Torkkuajan keston valinta asetuksista on selitetty osiossa "Hälytykset ja varoitukset" sivulla 69.

Kun napautat kellokuvaketta Tiedot ja kunto -alueella (⑦ Kuva 26 ja ⑧ Kuva 27), näkyviin avautuu näyttö, jossa näkyvät kaikki aktiivisen hälytyksen tiedot, ks. Kuva 37. Voit vierittää listaa ja napauttaa haluamaasi kohdetta laajentaaksesi sen ja lukeaksesi enemmän.



Kuva 37: Hälytysluettelo

Voit valita aktiivisen hälytyksen napauttamalla hälytyskohteen vieressä olevaa valintaruutua. Jos painat painiketta **Kuittaa**, hälytys yritetään poistaa luettelosta, jos hälytyksen pääsyy on poistettu. Kun kaikki hälytykset on poistettu, Tila-kuvakkeeksi palautuu oletusarvoinen vihreä sydän.

# <u>Virhe</u>

Toimintahäiriön ilmetessä ravistin antaa virheilmoituksen ja pysähtyy välittömästi, jotta näytteiden ja/tai yksikön vahingoittuminen voidaan välttää. Näyttö muuttuu kokonaan punaiseksi eikä käytön jatkaminen ole mahdollista. Näyttöön ilmestyy virheilmoitus virhekoodeineen, ks. Kuva 38.



Kuva 38: Virhenäyttö

Toimi ongelman suhteen seuraavasti ja yritä jatkaa käyttöä:

- 1. Merkitse muistiin virhenäytössä näkyvä virhekoodi.
- 2. Napauttamalla Kuittaa voit mykistää hälytysäänen.
- 3. Käynnistä ravistin uudelleen sammuttamalla se ensin ja käynnistämällä sitten uudelleen.
- 4. Jos virheilmoitus ei poistu, ota yhteys Thermo Fisher Scientificin asiakaspalveluun ja ilmoita virhenäytössä näkyvä virhekoodi.

# 3. 2. 3. Asetukset

Siirtymispalkin toinen kohde on Asetukset-kuvake. Alla oleva näyttö tulee näkyviin, kun napautat **Asetukset**. Asetukset-näytössä on enemmän painikkeita kuin näyttöikkunassa voi kerralla näyttää. Muut painikkeet tulevat näkyviin vierittämällä, ks. Kuva 39 oikea puoli.



Kuva 39: Asetukset-näyttö

# <u>Hälytykset ja varoitukset</u>

Hälytykset ja hälytysilmoitukset -näytössä voit määrittää, miten ja milloin hälytyksiä ja hälytysilmoituksia annetaan.



Kuva 40: Määritä Hälytykset ja hälytysilmoitukset -asetukset

Äänenvoimakkuus: Tällä säätimellä voit muuttaa hälytyksen tai hälytysilmoituksen yhteydessä kuuluvan äänen voimakkuutta. Vähennä äänenvoimakkuutta vetämällä liukusäädintä vasemmalle tai lisää sitä vetämällä liukusäädintä oikealle ja napauta sen jälkeen **Tallenna**. Tämä äänenvoimakkuusasetus vaikuttaa sekä hälytyksiin että hälytysilmoituksiin.



Kuva 41: Määritä kaiuttimen äänenvoimakkuus hälytyksille ja hälytysilmoituksille

Korkean lämpötilan hälytys / matalan lämpötilan hälytys (vain lämpötilaohjatut ravistimet): Näiden kahden asetuksen avulla voit määrittää lämpötilan asetusarvoon suhteutetun korkean ja matalan lämpötilan hälytyskynnyksen (ks. "Lämpötilan määritys" sivulla 64). Aina kun ravistimen kammion lämpötila laskee alle matalan lämpötilan kynnyksen tai ylittää korkean lämpötilan kynnyksen, ravistin antaa hälytyksen. Valitse lämpötilan kynnysarvo ja napauta Tallenna.



Kuva 42: Korkean ja matalan lämpötilan hälytyksen määritys

HUOMIO Aina, kun muutat lämpötilan asetusarvoa, hälytyskynnykset siirtyvät vastaavasti.

Korkean lämpötilan aiheuttama sammutus/ matalan lämpötilan aiheuttama sammutus (vain lämpötilaohjatut ravistimet): Näiden kahden asetuksen avulla voit määrittää lämpötilan kulloiseenkin asetusarvoon suhteutetun korkean ja matalan lämpötilan vaatiman laitteen sammutuksen turvasyistä (ks. "Lämpötilan määritys" sivulla 64). Kun ravistimen kammion lämpötila laskee alle matalan lämpötilan kynnyksen tai ylittää korkean lämpötilan kynnyksen, ravistin sammuu automaattisesti

kammiossa käsiteltävien näytteiden suojaamiseksi. Valitse kynnysarvo ja napauta **Tallenna**.



Kuva 43: Korkean ja matalan lämpötilan sammutuskynnykset

HUOMIO Aina, kun muutat lämpötilan asetusarvoa, virhekynnykset siirtyvät.

**Torkun aikakatkaisu**: Torkun aikakatkaisujakso määrittää, kuinka kauan hälytys pysyy mykistettynä, kun pääikkunan päällä näkyvän punaisen palkin torkkupainiketta painetaan (ks. kohtaa "Hälytys" sivulla 67). Voit asettaa torkkuajastimen 5, 10 tai 15 minuutille. Oletusarvo on 10 minuuttia. Napauta **Tallenna** tallentaaksesi muutokset.

Snooze Timeout	5min 🔨
	10min
	15min

Kuva 44: Hälytysten torkkuajastimen asetus

Poista kalibrointi-ilmoitukset käytöstä (vain lämpötilaohjatut ravistimet): Näiden asetusten avulla voit poistaa käytöstä ilmoitukset, jotka muistuttavat sinua säännöllisesti kalibroimaan ravistimen lämpötilamittauksen (ks. kohtaa "Lämpötilan kalibrointi" sivulla 126). Voit poistaa kalibrointimuistutukset käytöstä kokonaan tai vain siksi aikaa, kun ohjelma on käynnissä.

Disable Calibration Notifications Disable Calibration Notifications while a program is actively running Kuva 45: Poista kalibrointi-ilmoitukset käytöstä

### Pääsyn valvonta

Käytönvalvonta -näytössä voit määrittää järjestelmänvalvojan salasanan, jotta voit käsitellä prosesseja, joissa on rajoitetut käyttöoikeudet.

Ravistimen tila on oletusarvoisesti "Avoin tila", so. yksikön käyttämiseen ei tarvita salasanaa. "Suojattu tila" -tilaan siirtyminen edellyttää järjestelmänvalvojan salasanaa jokaiselta käyttäjältä, joka haluaa muuttaa ravistimen asetuksia. Kosketusnäytön oikeassa yläkulmassa näkyy tällöin Kirjaudu sisään -painike..

Seuraava vertailu tilojen "Avoin tila" ja "Turvatila" välillä osoittaa, milloin salasana tarvitaan.

Toiminto	Salasana tarvitaan avoimessa tilassa	Salasana tarvitaan turvatilassa
Nopeuden, ajan ja lämpötilan asetusarvojen määritys	Ei	Ei
Käynnistä ja pysäytä ravistin	Ei	Ei
Laske radat	Ei	Ei
Kuittaa hälytykset ja varoitukset	Ei	Ei
Katso tiedostoja ja tietonäyttöä	Ei	Ei
Näytä laitteen kunto	Ei	Ei
Näytä ja vie tapahtumaloki ja kaaviot	Ei	Ei
Torkkuhälytykset	Ei	Ei
Aja ohjelmia	Ei	Ei
Luo, muokkaa ja poista ohjelmia	Ei	Kyllä
Muuta näyttöasetuksia	Ei	Kyllä
Muuta valvonta-asetuksia	Ei	Kyllä
Muuta hälytys- ja hälytysilmoitusasetuksia	Ei	Kyllä
Muuta käytönvalvonta-asetuksia	Ei	Kyllä
Tuo tai vie ohjelmia	Ei	Kyllä
Palauta tehdasasetukset	Ei	Kyllä
Suorita kalibrointi	Ei	Kyllä
Asenna laiteohjelmistopäivitys	Kyllä	Kyllä

Taulukko 52: Salasanallisen kirjautumisen vaatimukset avoimessa ja turvatilassa

#### Graafinen käyttöliittymä

Kuva 46 näkyy Käytön valvonta -näyttö Avoin- ja Suojattu-tilassa.

< 余	Shake Lab 1 11:22:24AM Access Control		< *	Shake Lab 1 11:22:24AM Access Control	Logout Admin
¢.	<ul> <li>Open</li> <li>Secure Will require passcode</li> </ul>	access to	¢. (	Open Secure Will require passcode	access to
	change certain paran Change Admin Passcode	neters >		change certain param	neters
<b>.</b>	Save	0		Save	0
The start		0			U

Kuva 46: Pääsyn valvonta: Avoin tila (vasen) ja suojattu tila (oikea)

Kun napautat kohtaa **Suojattu** ja sen jälkeen **Tallenna**, sinua pyydetään antamaan järjestelmänvalvojan salasana, ennen kuin muutos voidaan vahvistaa. Kaikki yksiköt toimitetaan samalla tehtaalla esiasetetulla ja käsikirjaan painetulla järjestelmänvalvojan salasanalla.

Samoin käyttäjää, joka yrittää muuttaa salasanasuojattua asetusta, pyydetään antamaan järjestelmänvalvojan salasana.

HUOMIO Esiasetettu järjestelmänvalvojan salasana on "00000".

Vaihda järjestelmänvalvojan salasana -painikkeella voit muuttaa esiasetettua salasanaa. Jos napautat Vaihda järjestelmänvalvojan salasana -painiketta, sinua pyydetään antamaan ensin nykyinen salasana ja sen jälkeen uusi salasana ja vahvistamaan se vielä kerran. Uusi salasanasi on tallennettu -viesti ilmoittaa, että tallennus onnistui.

#### Kalibraatio

Kalibrointinäyttö (vain lämpötilaohjatut ravistimet) avaa joukon näyttöjä, joiden avulla voit kalibroida ravistimen tarkkaa lämpötilan säätöä varten. Kalibrointiprosessi on kuvattu kohdassa "Lämpötilan kalibrointi" sivulla 126.
### Hallintalaitteet

Säätimet-näytössä on asetukset, joiden avulla voit tehdä esiasetukset haluamillesi laitteen oletusarvoisille käyttöparametreille seuraavasti:

**Nopeuden asetuspiste:** Voit esiasettaa nopeuden asetusarvon välille 15–525 rpm. Valitse esiasetettu nopeus pyörittämällä valintakiekon valitsimia ja napauta **Tallenna**. Yksityiskohtaiset ohjeet Nopeuden asetusarvo -näytön käyttöön ovat osiossa "Määritä nopeus" sivulla 61 osiossa 3. 2. 1.

**Aikamoodi:** Voit käyttää ravistinta jatkuvassa tai ajastetussa tilassa. Jatkuvassa tilassa ravistin pysäytetään manuaalisesti käyttäjän harkinnan mukaan. Ajastetussa tilassa ravistimen käyttölaite pysähtyy automaattisesti, kun ajastimeen määritetty aika on kulunut loppuun. Ajastettua tilaa varten tilassa voit esiasettaa oletusarvoisen käyttöajan ja valita, näytetäänkö graafisessa käyttöliittymässä kulunut aika vai jäljellä oleva aika ravistimen ollessa käynnissä. Yksityiskohtaiset ohjeet Ajastettu tila -näytön käyttöön ovat osiossa "Määritä käyntiaika" sivulla 63 luvussa 3. 2. 1.

Lämpötilan asetusarvo (vain lämpötilaohjatut ravistimet): Tällä asetuksella voit määrittää oletuslämpötilan, jonka haluat näkyvän ravistimen näytössä käynnistyksen yhteydessä. Yksityiskohtaiset ohjeet Lämpötilan asetusarvo -näytön käyttöön ovat osiossa "Lämpötilan määritys" sivulla 64 luvussa 3. 2. 1.

Automaattinen uudelleenkäynnistys: Tämä toiminto käynnistää laitteen uudelleen normaalin käynnistä/pysäytä -käytön, ohjelman suorittamisen tai lämpötilan kalibrointitoiminnon aikana ilmenneen virtakatkoksen jälkeen. Jos automaattinen uudelleenkäynnistys on asetettu kohtaan **Ei**.

#### <u>Näyttö</u>

Näyttöasetukset antavat sinulle mahdollisuuden vaihtaa eri näyttövaihtoehtojen välillä. Kirkkaus: Säätääksesi näytön kirkkautta, käytä liukukytkintä tai +/- -painikkeita.



Kuva 47: Näytön kirkkauden säätö

Kieli: Vaihda näytön kieli napauttamalla Kieli -painiketta. Valitse haluamasi kieli valintakiekkoa kiertämällä ja napauta Tallenna.



Kuva 48: Valitse Näytön kieli

**Mittayksiköt (vain lämpötilaohjatut ravistimet):** Napauttamalla **Mittayksiköt** voit vaihtaa graafisen käyttöliittymän näytön kaikkien lämpötilalukemien lukemia Celsius °C -yksiköiden ja Fahrenheit °F -yksiköiden välillä.

Units of Measure	°C 🔨
	۴F

Kuva 49: Valitse lämpötilan näyttöyksikkö

**Päiväys**: Asettaaksesi päiväyksen napauta Päiväys -painiketta. Valitse haluamasi kuukausi-, päivä- ja vuosi valintakiekkoa kiertämällä ja napauta **Tallenna**.



Kuva 50: Määritä kuukausi, päivämäärä ja aika

Päiväyksen muoto: Asettaaksesi päiväysmuodon, napauta Päiväysmuoto -painiketta. Valitse haluamasi päivämäärän muoto (esimerkiksi KK/PP/VVV, näkyy muodossa 15.04.2015) ja napauta Tallenna.



Kuva 51: Määritä päivämäärämuoto

Aika: Asettaaksesi ajan ja aikaformaatin, napauta Aika -painiketta. Napauta **12 h** tai **24 h**, kierrä sen jälkeen valintakiekon tunti-, minuutti- ja AM/FM (vain 12 tunnin muoto) -valitsimia ja napauta **Tallenna**.



Kuva 52: Aseta aika ja valitse 12 tai 24 tunnin kello

Nukkumismoodi: Napauta Lepotila halutessasi määrittää ravistimen näytön siirtymään lepotilaan 15 minuutin käyttämättömyyden jälkeen. Lepotilassa näytössä näkyy Herätä napauttamalla -kehote.



Kuva 53: Ota lepotila käyttöön

**Alue**: Voit määrittää alueen, jolla laitetta käytetään, napauttamalla **Alue** -painiketta. Napauta Kaupunki, maa -tekstiruutua ja kirjoita kaupungin nimen ensimmäiset kirjaimet. Graafinen käyttöliittymä alkaa ehdottaa vastaavuuksia kolmen ensimmäisen merkin perusteella. Hyväksy jokin vastaavuus tai jatka koko nimen kirjoittamista ja napauta sen jälkeen **Tallenna** näyttönäppäimistöllä.

< *	Shake Lab 1 11:22.24AM Region										< ♠	Shake Lab 1 11:22:24AM Region	$\bigcirc$	
<b>≎</b> 。	Cit	y, Cour	ntry:	Enter the	city and d	country o	closest to	your tirr	e zone		⇔. (	City, Country:	Enter the city and country closest to your time zone New York, United States	
Q	W	E	R	т	Y	U	1	0	Ρ		Ê			
A	S	D	F	G	н	J	К	L	S	ave				
1	Z	X	С	V	В	N	М			1				
?!1:	23								-	<b>W</b>			Save	()

Kuva 54: Määritä alue

Laitteen nimi: Kun haluat antaa tai vaihtaa ravistimen nimen, napauta Yksikön nimi -painiketta. Napauta Yksikön nimi -tekstiruutua ja ala kirjoittaa. Kun olet valmis, valitse Tallenna näyttönäppäimistöllä.

< ♠	Sh 11 Un	ake La :22:24/ nit Nam	b 1 AM ie			2					< ♠	Shake Lab 1 11:22:24AM Unit Name	$\bigcirc$	
٥,	U	nit Na	me	Sha	ake La	ab 1				]	۵.			
Q	W	E	R	Т	Y	U	I.	0	Ρ			Unit Name:	Shake Lab 1	
А	S	D	F	G	Н	J	K	L	Sa	ive				
1	Z	X	С	۷	В	Ν	М	@	•	1	-		Save	0
?!1:	23							-	-	<b>.</b>			Save	0

Kuva 55: Anna ravistimelle nimi

Mukauta valikkoa: Voit mukauttaa siirtymispalkin kahta alakuvaketta napauttamalla Mukauta valikkoa -painiketta. Vedä haluamasi kuvake päänäytön alueelta korvattavaksi haluamasi kuvakkeen päälle. Vahvista napauttamalla **Tallenna**. HUOMIO Voit palauttaa siirtymispalkin tehdasasetukset milloin tahansa napauttamalla Nollaa valikko.



Kuva 56: Muokkaa siirtymispalkkia

#### Tiedostot ja info

Tässä näytössä näkyvät sarjanumero, graafisen käyttöliittymän sovellusohjelmisto ja sen versio, ravistimen pääohjaimen laiteohjelmisto ja sen versio, parametritiedosto ja sen versio sekä nykyinen käyttöjärjestelmä.

Voit myös palauttaa ravistelijan tehdasasetukset napauttamalla **Tehdasasetusten palautus**. Tehdasasetusten palautus edellyttää järjestelmänvalvojan salasanaa ja poistaa kaikki graafisen käyttöliittymän avulla tehdyt asetukset. Tehdasasetusten palautus ei poista tapahtumalokia.

< *	Shake Lab 1 11:22:24AM Files and Info		< *	Shake Lab 1 11:22:24AM Files and Info		
¢, ( ≝	Serial Number 123TSX990960 H.M.I. 4.1.0 Main Controller 7000 0612 V3 Parameters 50154439 V1	î	≎, ( ≝	Main Controller 7000 0612 V3 Parameters 50154439 V1 Board Support Package: 4.1.15-jethro-15.0		Î
Ē	Board Support Package: 4.1.15-jethro-15.0		Ē	Factory Reset	>	
<b>.</b>	Factory Reset	> 🗸		Firmware	>	Ļ
		()				()

Kuva 57: Tiedostot ja info

Kun vierität näyttöä alaspäin, näet **Laiteohjelmisto** -painikkeen Kuva 57 oikealla puolella. Voit asentaa ravistimeen uuden laiteohjelmiston napauttamalla tätä painiketta. Yksityiskohtaiset ohjeet uuden laiteohjelmiston asentamiseen on annettu kohdassa "4. 7. Laiteohjelmiston asennus" sivulla 130.

HUOMIO Uusi laiteohjelmisto tulee hankkia valtuutetulta huoltoteknikolta.

#### Kunnossapito

Huoltoasetukset ovat rajoitettuja ja niitä voivat käyttää valtuutetut huoltoteknikot.

## 3. 2. 4. Ohjelmat

Avaa ohjelmaluettelo tai luo uusi ohjelma napauttamalla siirtymispalkin **Ohjelmat**-kuvaketta. Voit luoda, muokata, poistaa, tuoda ja viedä ohjelman. Kuva 58 näkyy uuden laitteen (vasemmalla) ja jonkin aikaa käytössä olleen (oikealla) laitteen Ohjelmat-ikkuna. Käytössä oleva laite näyttää luettelon käyttäjän luomista ohjelmista.



Kuva 58: Ohjelmat

Jokaisessa ohjelmaluettelon kohdassa on silmäkuvake 💽, joka avaa Ohjelman pikakatselu -näytön, jossa on ohjelman tiedot, ja kynäkuvake, 🖍 jonka avulla ohjelma voidaan avata muokattavaksi.

#### Määritä kaikkien ohjelmien asetukset

Ohjelma-näytöstä pääset Lisäasetukset -painikkeella Lisäasetukset -näyttöön. Tässä näytössä voit määrittää, mitä tapahtuu, kun käynnistät ohjelman. Nämä asetukset koskevat kaikkia luomiasi ohjelmia.



Kuva 59: Määritä ohjelman asetukset

1. Napauta Lisäasetukset -painiketta.

- 2. Valitse **Näytä ajastin** -vaihtoehto, jos haluat määrittää, miten haluat ohjelman ajastimen näkyvän:
  - » Kulunut: kuinka kauan ravistin on ollut käynnissä Käynnistys-painikkeen painamisesta lähtien tai
  - » Jäljellä oleva: kuinka kauan ravistin pysyy käynnissä, kun ajastimen aika on kulunut loppuun
- Jos haluat, että ravistin käynnistyy heti ohjelman käynnistämisen jälkeen, ota Käynnistä ravistin ohjelman latauksen yhteydessä -asetus käyttöön. Oletusmenettely on, että ohjelman suorittaminen edellyttää Käynnistä -painikkeen napautusta.

HUOMIO Käynnistä **ravistin ohjelman latauksen yhteydessä** toimii odotetulla tavalla vain, jos ravistimen kansi on kiinni. Kun kansi on suljettu, toiminta käynnistyy automaattisesti.

4. Napauta Tallenna.

#### Ohjelman luonti

Voit luoda ja tallentaa jopa 99 ohjelmaa.

- 1. Napauta Luo uusi-painiketta.
- 2. Anna ohjelman nimi.





3. Syötä ohjelma lämpötila (vain lämpötilaohjatut ravistimet), nopeus ja käyntiaika tunteina ja minuutteina.

Jos haluat poistaa lämpötilaohjatun ravistimen lämpötilan säädön käytöstä, jätä Lämpötila-kenttä tyhjäksi.

- Kun haluat lisätä vaiheen ohjelmaasi, vieritä näyttöä alaspäin ja napauta Lisää vaihe -painiketta.
- 5. Napauta **Tallenna** tallentaaksesi ohjelman.

#### Ohjelman muokkaus

- 1. Naputa lyijykynäkuvaketta sen ohjelman 🖍 vieressä, jota haluat muokata.
- 2. Muokkaa halutut kentät. Napauta **Tallenna**. Ohjelma tallennetaan uusien muutoksien kanssa.
- 3. Voit lisätä vaiheen valitsemalla Lisää vaihe näytön alaosassa. Vieritä näyttöä nähdäksesi Lisää vaihe -painikkeen, jos vaiheita on enemmän kuin 3.

HUOMIO Jos sekoittamisen, jossa on useampia vaiheita, hälytys tai virheviesti tapahtuu, ravistin pysähtyy automaattisesti. Jos varoitus annetaan, sekoitus jatkuu.



Kuva 61: Lisää vaihe

#### Ohjelman poistaminen

- 1. Naputa lyijykynäkuvaketta sen ohjelman vieressä, jota haluat muokata,
- 2. Napauta **Poista**. Ikkuna ilmestyy ja kysyy varmennusta.



Kuva 62: Poista ohjelma

3. Napauta Poista. Valittu ohjelma poistetaan.

#### Käytä ohjelmaa

- 1. Käyttääksesi olemassa olevaa ohjelmaa, valitse ohjelma jota haluat käyttää.
- Jos haluat nähdä pikakatsauksen ohjelmasta, napauta silmäkuvaketta ohjelman vieressä.
- 3. Pikakatsausikkuna ilmestyy.



Kuva 63: Ohjelman pikakatsaus

4. Napauta Käytä ladataksesi ohjelman Kotinäytölle.

Valituista **Automaattinen toiminta** -asetuksista riippuen ohjelma käynnistyy välittömästi tai se käynnistyy sitten, kun **Käynnistys** -painiketta napautetaan kotinäytöllä.

5. Kun ohjelma on valmis, seuraava ikkuna ilmestyy. Napauta OK.



Kuva 64: Ohjelma valmis

#### Ohjelmien tuonti

Voit tuoda yhdellä ravistimella luotuja ohjelmia toiseen ravistimeen. Tämä edellyttää USB-aseman liittämistä ravistimen USB-porttiin.

Seuraavassa näytössä näkyvät kaikki USB-asemassa tunnistetut ohjelmat.



Kuva 65: Vallitse tuotavat ohjelmat

Valitse ohjelmat, jotka haluat tuoda. Napauta Tuo.

Kun ohjelmat on onnistuneesti tuotu, seuraava ikkuna ilmestyy. Napauta **OK**. Voit nyt postaa USB:n.



Kuva 66: Ohjelman tuonti valmis.

#### Ohjelmien vienti

Voit viedä ohjelmia yhdestä ravistimesta toiseen. Varmista, että USB-asema on kytketty.

Seuraava näyttö näkyy jos USB on yhdistetty:



Kuva 67: Valitse ohjelmat vientiä varten

Valitse ohjelmat, jotka haluat viedä. Napauta Vie.

Seuraava ikkuna ilmestyy kun vienti on valmis. Napauta OK. Voit nyt postaa USB:n.



Kuva 68: Ohjelman vienti valmis

### <u>Tapahtumaloki</u>

Kolmas välilehti navigaatiopaneelissa on tapahtumaloki, joka sisältää talletetut käyttäjä- ja järjestelmätapahtumat. Tapahtumaloki-näyttö avautuu napauttamalla valikkorivin Tapahtumaloki-kuvaketta, ks. Kuva 69.



Kuva 69: Tapahtumaloki

Näyttö näyttää 50 viimeisintä tapahtumaa, kunkin päivämäärä- ja aikaleimoineen.

Lisätietoja yksittäisestä tapahtumasta voi katsoa valitsemalla tapahtuman luettelosta.

Tapahtumat voidaan suodattaa tyypin mukaan seuraaviin luokkiin:

- hälytykset
- varoitukset
- asetusten muutokset
- käynnistys/pysäytystoiminnot
- ohjelman käytöt
- kannen avaustoiminnot (vain lämpötilaohjatut ravistimet)

Kun suodatin valitaan, oikean puolen Näytä kaikki -painike vaihtuu Suodatin päällä -painikkeeksi, ks. Kuva 70.

	Shake Lab 1 11:22:24AM Event Log	Ø	)		
	Date/Time	Event	Filter ON	$\mathbf{v}$	
٥.	5/12/15 11:34:06AM	Setpoint Change		$\mathbf{\sim}$	
	5/12/15 11:34:06AM	High Temp Alarm		$\boldsymbol{\sim}$	
	5/12/15 11:34:06AM	Unit Stopped		$\sim$	
Ļ	5/12/15 11-24-06AM	Unit Started		$\checkmark$	$\sim$
	Export	Log	View Cha	rt	í

Kuva 70: Suodatettu tapahtumaloki

#### Tapahtumalokin vienti

1. Valitse vietävät tapahtumat avattavasta Vietävät tapahtumat -luettelosta. Valitse lokin tai raportin vientimuoto.



Kuva 71: Valitse tapahtumaloki vietäväksi

2. Ennalta määritetty tai kustomoitu päiväysalue voidaan valita.

HUOMIO Voit viedä tapahtumia vain kuuden viime kuukauden ajalta.

 Vienti voidaan tehdä käyttäen USB-muistia. USB-muistitikku täytyy olla kytkettynä laitteeseen lokin tai raportin tallentamiseksi. Napauta Vie loki -painiketta ladataksesi lokin tai raportin.



Kuva 72: Aseta USB-muisti vientiä varten.

## <u>Kaaviot</u>

Kaavioissa näkyvät ajan funktioina nopeus- tai (vain lämpötilaohjatut ravistimet) lämpötilatietojen käyrät. X-akselilla näkyy aika ja Y-akselilla nopeus tai lämpötila.

Lämpötilakaavion avulla voit piirtää asetusarvon, todellisen lämpötilan ja ympäristön lämpötilan ajan funktiona. **Muokkaa**-painikkeella voit valita, mitkä kolme käyrää lämpötilakaavioon sisällytetään. Lisäksi katselualueen voi muuttaa 1 päiväksi, 7 päiväksi tai käyttäjän määrittämäksi määräksi päiviä joko täyden ajalle 24 tunnin ajan tai valituille tunneille kunakin päivänä.

**Päivitä**-painike päivittää kaavionäytön asetustesi mukaiseksi ja näyttää kaikki uudet tiedot, jotka on tallennettu sen jälkeen, kun kaavio viimeksi ladattiin tai päivitettiin.

**HUOMAUTUS** Lämpötilahälytyksen sattuessa graafisessa käyttöliittymässä näkyvä lämpötila vaihtuu välittömästi ravistimen kammion todelliseen ilman lämpötilaan, jotta näytteitä voitaisiin suojata mahdollisimman hyvin. Tällöin kaavio-ominaisuudessa esitetyssä lämpötilassa voi näkyä hyppy.



Kuva 73: Nopeus- ja lämpötilakaaviot

Voit asettaa USB-aseman ravistimen USB-porttiin ja ladata nopeus- tai lämpötilakaavion tiedot napauttamalla **Vie karttatiedot**.

#### Kaavion muokkaus

1. Aloita kartan muokkaaminen napauttamalla Muokkaa.



Kuva 74: Kaavion muokkaus Valitse tallennettavaksi lämpötila tai nopeus

2. Valitse näytettävä kaavio napauttamalla Lämpötila tai Nopeus.

- 3. Vieritä näyttöä alaspäin niin, että näyttöön tulevat Päivämääräväli -asetukset.
- 4. Valitse napauttamalla 1 päivä, 7 päivää tai **Mukautettu** mukautetulle päivämäärävälille.
- 5. Valitse aloituspäivämäärä pyörittämällä valintakiekon valitsimia ja napauta **Aseta** aloituspäivä -painiketta.
- Valitse lopetuspäivä pyörittämällä valintakiekon valitsimia ja napauta Aseta lopetuspäivä -painiketta.



Kuva 75: Valitse kaavion päivämääräväli

- 7. Vieritä näyttöä alaspäin niin, että näyttöön tulevat Aikaväli -asetukset.
- 8. Valitse 24 tunnin päivä napauttamalla tai napauta **Mukautettu**, jos haluat kaavioon tietyn ajanjakson ja
  - a. Valitse päivittäinen aloitusaika pyörittämällä valintakiekon valitsimia ja napauta **Aseta aloitusaika** -painiketta.
  - b. Valitse päivittäinen lopetusaika pyörittämällä valintakiekon valitsimia ja napauta Aseta lopetusaika -painiketta.

**HUOMIO** Valitsemalla aloitus- ja lopetusajat voit valita kaavioon tarkan kahden ajankohdan välisen ajanjakson. Jos käytät 24 tunnin päivä -asetusta, kaavion aloitusajaksi tulee klo 00:00 ja päättymisajaksi 23:59 joka päivälle.



Kuva 76: Valitse kaavioon tulevat päivän tunnit

9. Napauta Tallenna tallentaaksesi muutokset tai räätälöidyt asetukset.

Tarvikkeet

## 3.3. Tarvikkeet



Fyysinen ja biologinen haitta johtuen rikkinäisistä astioista. Väärin asennetut apuvälineet voivat johtaa rikkinäiseen lasiin ja läikkyneisiin näytteisiin.



Varmista, että apuvälineet on oikein asennettu käyttäen oikeita työkaluja ja ruuveja.

Pidä huolta, että apuvälineet sopivat alustaan mielekkäästi. Käytä aina sopivan kokoista astiaa ja lisävarustetta.



Viiltohaavat terävistä reunoista.

NOUDATA VAROVAI-SUUTTA Ole varovainen, kun käsittelet alustoja ja muita apuvälineitä.



NOUDATA VAROVAI-SUUTTA Sormet voivat jäädä puristuksiin liikkuvaan alustaan. Älä koskaan laita sormiasi liikkuvan alustan päälle tai alle.

**HUOMIO** On sinun vastuullasi varmistaa, että apuvälineet on asennettu asianmukaisesti.

Asennustoimenpiteet, jotka kuvataan, koskevat kaikkia ravistinmalleja, jotka on listattu tässä käyttöoppaassa, jos muuta ei mainita.

Käytä aina oikeita työkaluja, erityisesti tarvikkeen mukana toimitettuja työkaluja. Jos kadotat joitakin näistä esineistä, voit tilata luettelossa mainittuja varaosasarjoja ("Tarvikkeet" sivulla 17). Älä käytä muita kuin määritettyjä työkaluja tai muita kuin toimitukseen sisältyviä ruuveja.

## 3. 3. 1. Alustan asennus

NOUDATA VAROVAISUUTTA Täydellinen luettelo kunkin ravistimen kanssa yhteensopivista alustoista on osiossa "1. 2. 1. Alustat" sivulla 18. Lämpötilaohjattuja ravistinmalleja saa käyttää vain tehtaalla asennettujen alustojen kanssa.

**NOUDATA VAROVAISUUTTA** Viiltohaavat terävistä reunoista. Kun poistat alustaa, ota kiinni alustasta sen alapuolelta. Älä nosta alustaa puristimista, jotka on siihen kiinnitetty. Ole varovainen, kun käsittelet alustoja ja muita apuvälineitä.

NOUDATA VAROVAISUUTTA Solaris 4000 I / 4000 R -ravistimen liikkuva 11x14-alusta voi jättää kädet tai sormet puristuksiin. Älä käytä 11x14 -alustaa Solaris 4000 I / 4000 R -ravistimessa.

NOUDATA VAROVAISUUTTA Käytä vain ruuveja ja lukitustyökalua, jotka tulevat mukana. Toisten ruuvien tai väärien lukitustyökalujen käyttö johtaa väärään asennukseen ja voi mahdollisesti vahingoittaa ravistinta ja apuvälineitä.

Asenna aina alusta **kaikilla** ruuveilla.

Käytä aina oikeata lukitustyökalua:

- 3/16 tuuman lukitustyökalu (GT530066) kaikille Solaris 2000 ja Solaris 2000 I/R - ja 4000 I/R -ravistimien alustoille
- 7/32 tuuman lukitustyökalu (GT530080) kaikille Solaris 4000 -ravistimenalustoille

## <u>Yleisalustat</u>

Jokaisessa Solaris 2000- ja 4000-ravistimessa on yksi yleisalusta, ruuvit ja lukitustyökalu. Voit ostaa omiin käyttötarkoituksiisi sopivia lisäalustoja. Täydellinen luettelo kunkin ravistimen kanssa yhteensopivista alustoista on osiossa "1. 2. 1. Alustat" sivulla 18.

2000 I -, 2000 R -, 4000 I - ja 4000 R -ravistimissa on kussakin ravistimeen valmiiksi asennettu alusta. Alustaa tai ruuveja ei ole saatavana erillisinä osina.



Kuva 78: Yleisalusta asennettuna Solaris 2000 (vasemmalla) ja Solaris 4000 (oikealla) -ravistimeen

1. Aseta alusta varovasti vaakatasossa ravistimen ja sen kiinnityspisteiden yläpuolelle.

Solaris 2000 -ravistimessa on 3 kiinnitysreikää. Solaris 4000 -ravistimessa on 4 kiinnitysreikää.

Solaris 2000 I/R - ja 4000 I/R -ravistimissa on 3 kiinnitysreikää.

2. Kiristä ruuvit varoen kiinnittääksesi alustan ravistimeen. Kun lukitustyökalu alkaa taipua, lopeta kiristäminen.

### Dual Stack-alustat

HUOMIO Kaksitasoiset alustat eivät ole yhteensopivia Solaris 2000 I/R - ja 4000 I/R -ravistimen kanssa.



Kuva 79: Dual Stack -alustan kokoaminen

Yhdistä ylempi ja alempi alusta kiinnittämällä ne 4 pilariin joka kulmassa. Kiinnitä pilarit sopivilla ruuveilla ylä- ja ala-alustoihin.

Kiristä ruuvit varoen yhdistääksesi pilarin ja alustat. Kun lukitustyökalu alkaa taipua, lopeta kiristäminen.

Ala-alustassa on kiinnitysreiät ravistimeen liittämiseen. Huolehdi, että ala-alusta on pohjapuolella alustaa asentaessasi.

#### Tarvikkeet

HUOMIO Ennen kuin asetat astioita alustarakenteelle, tee viimeinen tarkistus varmistaaksesi, ettei alustarakenne huoju.



Kuva 80: Dual stack -alusta Solaris 4000 -ravistimella

1. Aseta varovasti koottu dual stack-alusta vaakasuunnassa ravistimen ja sen alustan kiinnityslevyn yläpuolelle.

Alustoissa Solaris 2000:lle on 3 kiinnitysreikää, alustoissa Solaris 4000:lle on 4 kiinnitysreikää.

2. Kiristä ruuvit varoen kiinnittääksesi alustan ravistimeen. Kun lukitustyökalu alkaa taipua, lopeta kiristäminen.

### 3. 3. 2. Koepullopidikkeen ja astian asennus

NOUDATA VAROVAISUUTTA Rikkinäisistä tai vuotavista astioista johtuva biologinen haitta. Väärin asennetut apuvälineet voivat johtaa läikkyneisiin näytteisiin. Varmista, että apuvälineet on oikein asennettu käyttäen oikeita työkaluja ja ruuveja. Asennukseen tarvitaan ruuvitaltta (PH 2), jonka varren pituus on 150 mm / 6 tuumaa. (Tuotenumero 75004131). Pidä huolta, että apuvälineet sopivat alustaan mielekkäästi. Käytä aina sopivan kokoista astiaa ja lisävarustetta. Astioiden täytyy olla virheettömiä ja oikein asennettu.

NOUDATA VAROVAISUUTTA Viiltohaavat terävistä reunoista. Ole varovainen, kun käsittelet alustoja ja muita apuvälineitä.

NOUDATA VAROVAISUUTTA Pidike voidaan kiinnittää turvallisesti ainoastaan toimitukseen sisältyvien ruuvien avulla. Asenna toimitukseen sisältyvät ruuvit pidikkeen kaikkiin tähän tarkoitukseen varattuihin reikiin.

HUOMIO Jos ruuvit ovat liian pitkät, vaihdettavan alustan vakaus ja toiminta saattavat heikentyä. Varmista, että kiinnitykseen käytetään ainoastaan toimitukseen sisältyviä ruuveja.

## Puristimet

Jokainen puristin koostuu astiaklipsistä, yhdestä tai kahdesta jousesta riippuen puristimen/pullon koosta, ja ruuveista alustaan kiinnitykseen. Käytä vain puristimen mukana tulevia ruuveja.



1) Jousiputki 2) Jousi 3) Ruuvit 4) Klipsin jalka 5) Jousi

Kuva 81: Puristimen yksityiskohdat 2 jousella

- 1. Jos on tarpeen, kiinnitä jousi klipsin jalkoihin kuten kuvissa.
- Kumiset jousiputket sijaitsevat klipsin jalkojen välissä, kuten kuvassa. Joissakin pidikkeissä on kaksi jousta. Toinen jousi asennetaan klipsirakenteen pohjan ympäri kun se on kiinnitetty alustaan.
- 3. Kiinnitä klipsirakenne alustaan mukana tulevilla ruuveilla.

## <u>Astia</u>

 Aseta varovasti haluttu astia puristimeen ensin vetäen puristimen jousen niin pitkälle, että astian pohja menee hyvin puristimeen. Liu'uta varovasti astia oikeaan asentoon ja, kiinnitä se hyvin puristimen leveämpään pohjaan. Jousi pitää astian kaulan tiukasti paikoillaan.



2. Varmista, että kaikki astiat ovat hyvin kiinni, ennen kuin panet virran päälle.

Astiassa tulee mahdollisuuksien mukaan sulkea tulpalla vaarallisten aineiden läikkymisen estämiseksi sekoituksen aikana.

## 3. 3. 3. Nelikulmaisen pidikkeen asennus

NOUDATA VAROVAISUUTTA Rikkinäisistä tai vuotavista astioista johtuva biologinen haitta. Väärin asennetut apuvälineet voivat johtaa läikkyneisiin näytteisiin. Varmista, että apuvälineet on oikein asennettu käyttäen oikeita työkaluja ja ruuveja. Asennukseen tarvitaan ruuvitaltta (PH 2), jonka varren pituus on 150 mm / 6 tuumaa. (Tuotenumero 75004131). Pidä huolta, että apuvälineet sopivat alustaan mielekkäästi. Käytä aina sopivan kokoista astiaa ja lisävarustetta. Astioiden täytyy olla virheettömiä ja oikein asennettu.

NOUDATA VAROVAISUUTTA Viiltohaavat terävistä reunoista. Ole varovainen, kun käsittelet alustoja ja muita apuvälineitä.

## Nelikulmaisen pidikkeen tiedot

Jokaisessa nelikulmaisessa pidikkeessä on pohjapaneeli, jossa on kiinnitysreiät, ja neljä jousikuormitettua pidikejalkaa. Toimitukseen sisältyvät myös ruuvit, joilla pidike kiinnitetään alustaan. Käytä vain puristimen mukana tulevia ruuveja.



1) Kiinnitysreiät 2) Jousi 3) Pidikejalka 4) Ruuvit

Kuva 82: Nelikulmaisen pidikkeen tiedot

- 1. Kiinnitä tarvittaessa kaikki jouset 2 pidikejalkoihin 3, ks. Kuva 82.
- 2. Aseta pidike alustalle ja kohdista kiinnitysreiät ①.
- 3. Kiinnitä pidikerakenne alustaan toimitukseen sisältyvillä ruuveilla ④.
- 4. Työnnä nelikulmainen astia paikalleen ylhäältä.
- 5. Varmista, että astia on pohjapaneelin päällä ja että se on tukevasti pidikkeessään.

## 3. 3. 4. Koeputkitelineen asennus

NOUDATA VAROVAISUUTTA Viiltohaavat terävistä reunoista. Ole varovainen, kun käsittelet alustoja ja muita apuvälineitä.

Jokainen koeputkiteline koostuu kahdesta lukitusnupillisesta (metallilankakehyksisestä tai peltisestä) asennuskiinnikkeestä ja kunkin koeputkitelineen kahdesta siipiruuvista, joilla teline kiinnitetään alustaan. Jokainen koeputkiteline osineen toimitetaan esikoottuna, ja toimitukseen sisältyvät koeputkiteline ja vaahtomuovityyny tai -tyynyt. Käytä aina telineen mukana toimitettuja ruuveja.



Kuva 83: Koeputkiteline osineen, mukana metallilankaiset asennuskiinnikkeet



Kuva 84: Koeputkiteline osineen, mukana peltiset asennuskiinnikkeet

- 1. Kiinnitä asennuskiinnikkeet alustaan siipiruuveilla. Kiristä sormiruuveja, kunnes ne ovat käsivaralta tiukat.
- 2. Työnnä koeputkiteline asennuskiinnikkeisiin.
- 3. Aseta sopivaan kulmaan käyttäen lukitusnuppeja. Kiristä lukitusnuppeja, kunnes ne ovat käsivaralta tiukat.

# 3. 3. 5. Kallistuskulmaltaan säädettävän koeputkitelineen pidikkeen asennus

NOUDATA VAROVAISUUTTA Viiltohaavat terävistä reunoista. Ole varovainen, kun käsittelet alustoja ja muita apuvälineitä.

Koeputkipidikkeitä ja koeputkitelineitä on saatavana erikokoisina, ks. luku "1. 2. 5. Kallistuskulmaltaan säädettävien koeputkitelineiden pidikkeet" sivulla 28. Kaikki koeputkitelinepidikkeet voidaan säätää seitsemään asentoon, jolloin ne kääntyvät asennosta 0° kumpaankin suuntaan ja voidaan lukita asentoihin 15°, 30° ja 45°.

Käytä aina telineen pidikkeen mukana toimitettuja ruuveja.



Kuva 85: Kallistuskulmaltaan säädettävä koeputkitelinepidike ja koeputkiteline asennettuina

- 1. Kiinnitä telineen pidikkeet ① ravistimeen toimitukseen sisältyvillä ruuveilla.
- 2. Levitä metallikielekkeitä ③ telineen pidikkeen kääntöalustan kummassakin päässä ⑥ ja työnnä koeputkiteline ② paikalleen.
- 3. Kiinnitä koeputkiteline 2 vapauttamalla varovasti kielekkeet 3.
- 4. Vedä pidikkeen kummankin pään lukitustappien nuppeja 6 ulospäin. Tappien lukitus avataan kääntämällä nuppia 1/4 kierrosta.
- Käännä telineen pidikkeen ① kääntöalusta ⑥ haluamaasi kulmaan 15°, 30° tai 45°.
- Asenna lukitustappi (5) takaisin kohdistusreikään (4) ja lukitse tappi kääntämällä nuppia 1/4-kierrosta.
- 7. Kun haluat irrottaa telineen, levitä metallikielekkeitä ③ kääntöalustan ⑥ kummassakin päässä ja nosta koeputkiteline ① pois.

## 3. 3. 6. Mikrolevyn / Deepwell-levyn pidikkeen asennus

NOUDATA VAROVAISUUTTA Viiltohaavat terävistä reunoista. Ole varovainen, kun käsittelet alustoja ja muita apuvälineitä.



- 1) Deepwell-tarjotin
- 2 Ruuvi
- ③ Mikrotarjottimen puristin

Kuva 86: Mikrolevy / Deepwell-levy osineen

- 1. Aseta mikrotarjottimen kehys alustaan.
- 2. Kiinnitä mikrotarjottimen kehys alustaan käyttäen ruuveja, jotka tulevat mukana mikrotarjotin-/deepwell-tarjotin setin mukana.
- 3. Aseta mikrotarjotin tai deepwell-tarjotin mikrotarjottimen kehykseen.
- 4. Varmista, että mikrotarjotin tai deepwell-tarjotin istuu tiukasti nostamalla sitä kevyesti.

## 3. 3. 7. Säädettävän astianpidikkeen asennus

NOUDATA VAROVAISUUTTA Viiltohaavat terävistä reunoista. Ole varovainen, kun käsittelet alustoja ja muita apuvälineitä.





① Kiinnitysreiät ② Säätötangot ③ Lukitusnupit ④ Alustan sivulevy ⑤ Liukumaton matto
 ⑥ Ruuvit

Kuva 87: Yksittäisen säädettävän astianpidikkeen asennus

Tarvikkeet





① Kiinnitysreiät ② Säätötangot ③ Lukitusnupit ④ Alustan sivulevy ⑤ Liukumaton matto
 ⑥ Ruuvit

Kuva 88: Kahdelle korkeudelle säädettävä astianpidike osineen

Suorita asennus seuraavasti:

- 1. Nosta liukumatonta mattoa niin, että kiinnitysreiät tulevat näkyviin (kohteet Kuva 88).
- 2. Aseta säädettävä astianpidike alustalle ja kohdista kiinnitysreiät.
- 3. Kiinnitä säädettävä astianpidike alustaan kiinnikkeen mukana (© toimitetuilla ruuveilla.
- 4. Aseta liukumaton matto (5) säädettävän astianpidikkeen pohjalle.
- 5. Aseta astia(t) liukumattomalle matolle .
- 6. Löysää lukitusnupit 3.
- 7. Säädä säätötankojen asentoa 2 siten, että astiat ovat kunnolla kiinni.
- 8. Kiristä lukitusnupit 3.
- 9. Varmista, että astia on kunnolla kiinnitetty nostamalla sitä varovasti.

#### 100 / 140

## 3. 3. 8. Dekantterilasitelineen asennus

NOUDATA VAROVAISUUTTA Viiltohaavat terävistä reunoista. Ole varovainen, kun käsittelet alustoja ja muita apuvälineitä.



Kuva 89: Dekantterilasiteline osineen

- 1. Aseta pohjapaneeli 2 alustalle.
- 2. Aseta dekantterilasiteline ① pohjapaneelin päälle.
- 3. Kohdista kiinnitysreiät alustaan.
- 4. Kiinnitä dekantterilasiteline tuotteen mukana toimitetuilla ruuveilla 3.

## 3. 3. 9. Erotussuppilon pidikkeen asennus

NOUDATA VAROVAISUUTTA Viiltohaavat terävistä reunoista. Ole varovainen, kun käsittelet alustoja ja muita apuvälineitä.

HUOMIO Varmista, että asennuksessa käytetään oikeita ruuveja. Toimitukseen sisältyvät kahdet erilaiset ruuvit. Pystytuen kiinnittämiseen tulee käyttää pidempiä ruuveja, joissa on punainen kierrelukite. Lyhyempiä ruuveja, joissa ei ole kierrelukitetta, saa käyttää vain erotussuppilon pidikkeen kiinnittämiseen alustaan.

Erotussuppilon pidike voidaan asentaa ravistimen alustaan kahteen perusasentoon: vaakasuoraan ja pystysuoraan.



Kuva 90: Erotussuppilon pidikkeen kiinnitysasennot

Kuva 90 vaakasuoraa asentoa ③ voidaan säätää korkeussuunnassa siten, että saavutetaan kohdan ④ kallistettu asento. Lisäksi erotussuppilon pidike voidaan yhdistää pystytelineeseen (saatavana lisätarvikkeena), jolloin tuloksena ovat pystysuorat tyypit ① ja ②. Lisäksi pidike voidaan asentaa siten, että suppilon hana osoittaa alaspäin ① tai päinvastoin, ks. ylösalaisin olevaa järjestelyä kohdassa ②.

## Pystyasennuksen valmistelu

Kuva 91 näkyvät erotussuppilon pidike 🛞 ja lisätarvikkeena pystyjalusta 🕲 vierekkäin.



Kuva 91: Erotussuppilon pidike (vasemmalla) ja lisätarvikkeena toimitettava pystyjalusta (oikealla)

- 1. Aseta erotussuppilon pidike pystysuuntaisen kiinnikkeen pitkään päähän, ks. kohdat ① ja @ Kuva 90.
- 2. Kohdista seuraavaksi neljä pystyjalustan suppilon kiinnitysreikää (kohta @ Kuva 91):
  - a. Käytä vakiomallisen pystyasennuksen kiinnitysreikiä (3 Kuva 91) halutessasi asentaa suppilon hanan osoittamaan alaspäin.
  - b. Käytä ylösalaisin tulevan pystyasennuksen kiinnitysreikiä (④ Kuva 91) halutessasi asentaa suppilon hanan osoittamaan ylöspäin.
  - Kiinnitä erotussuppilon pidike pystyasentoon toimitukseen sisältyvillä ruuveilla (6 Kuva 91).

#### Asennus alustalle

- 1. Aseta erotussuppilon pidike alustalle.
- 2. Kohdista kiinnitysreiät alustaan.
- 3. Kiinnitä erotussuppilon pidike alustaan erotussuppilon pidikkeen mukana toimitetuilla ruuveilla (6 Kuva 91 ja Kuva 92).

#### Suppilon asennus



 Erotussuppilo
 Suppilon takapidike
 Suppilon etupidike
 Lukitusnuppi, korkeussäätö
 Lukitusnuppi, pituussäätö
 Ruuvit ilman kierrelukitetta

Kuva 92: Erotussuppilon pidike osineen

- 1. Löysää pituussäädön lukitusnuppia (5) ja liu'uta suppilon etupidike (3) kokonaan etuasentoon.
- 2. Aseta suppilon takaosa ① suppilon takapidikkeeseen ②.
- 3. Liu'uta suppilon etupidikettä ③ taaksepäin ja aseta suppilon varsi suppilon etupidikkeen aukkoon ③.
- Säädä suppilon etupidike ③ sopivalle korkeudelle ja kiinnitä se kiristämällä korkeussäädön lukitusnuppi ④. Voit kallistaa erotussuppilon vinoon korkeussäätöä muuttamalla.
- 5. Kiinnitä suppilon etupidike 3 kiristämällä pituussäädön lukitusnuppia 5.

## 3. 3. 10. Tarviketason asennus

NOUDATA VAROVAISUUTTA Viiltohaavat terävistä reunoista. Ole varovainen, kun käsittelet alustoja ja muita apuvälineitä.



Kuva 93: Tarviketaso osineen

Suorita asennus seuraavasti:

- 1. Aseta tarviketaso 2 alustalle.
- 2. Kohdista kiinnitysreiät alustaan.
- 3. Kiinnitä tarviketaso 2 ruuveilla 3, jotka on toimitettu tarviketason.
- 4. Aseta kumimatto 1 tasolle 2.
- Aseta pullot tasolle ja kiinnitä ne toimitukseen sisältyvillä kumirenkailla @, ks. Kuva 94.

HUOMIO Kumirenkaiden käyttöä suositellaan yli 100 rpm:n nopeuksille.



Kuva 94: Tarviketaso ja asennetut säiliöt ja kumirenkaat

HUOMIO Varo venyttämästä kumirenkaita liikaa kiinnittäessäsi säiliöitä. Kun venytät kutakin asennettavaa kumirengasta, huomaat, että joustavuus loppuu tietyssä pisteessä, minkä jälkeen kumirenkaat eivät enää veny. Kun olet saavuttanut tämän pisteen, älä yritä venyttää kumirenkaita enempää, koska muutoin rengas katkeaa.

## 3. 3. 11. Liimakiinnitteisen maton asennus ja käyttö

Liimakiinnitteiset matot soveltuvat leveisiin, tasaisille säiliöille, kuten erlenmeyerpulloille, fernbach-pulloille, reagenssipulloille, mikrolevyille, petrimaljoille, soluviljelypulloille, mittapulloille ja dekantterilaseille. Yleensä voidaan käyttää kaikkia astioita, jotka pysyvät pystyssä, mutta kun niiden pohjapinta-alan ja korkeuden suhde pienenee, niin pienenee myös niiden suurin mahdollinen ravistelunopeus. Korkeat säiliöt, joilla on pieni pohjapinta-ala, eivät ole yhtä sopivia, ks. Kuva 95. Suuret erlenmeyerpullot (esim. 3000 ml) kestävät paremmin kuin pienet (esim. 100 ml).



Kuva 95: Liimakiinnitteinen matto: Sopivat pullomuodot

Saatavana on kahdenlaisia liimakiinnitteisiä mattoja:

- Kevyesti kiinnittyvä matto (läpinäkyvä) 75004111 ja 75004117
  - » helppo käsitellä, koska tahmeus on vähäinen
  - » sopii erityisesti dekantterilaseille
  - » sopii kaikille astioille korkeintaan nopeudella 200 rpm (kudosviljelykäyttö)
- Tiukasti kiinnittyvä matto (musta läpinäkyvä) 75004126 ja 75004127
  - » luotettava astian kiinnitys, koska tahmeus on suuri
  - » sopii kaikille astioille
  - » enimmäisnopeus 400 rpm (enintään 500 ml:n erlenmeyerpullo)

## <u>Työkalut</u>

Kiinnitystela, joka näkyy Kuva 96, toimitetaan liimakiinnitteisten mattojen lisävarusteena. Se on tarkoitettu apuvälineeksi, jonka avulla voit painaa liimakiinnitteisen maton kiinni alustaan.



Kuva 96: Liimakiinnitteisen maton kiinnitystela

#### Kuivamaton asennus

- 1. Varmista, että ravistimen alusta on puhdas ja kuiva.
- 2. Irrota suojakalvo liimamaton toiselta puolelta.
- 3. Kiinnitä matto liimapuoli tukevasti alustaan.
- 4. Poista päällimmäinen suojakalvo.

#### Märkämaton asennus

- 1. Irrota suojakalvo liimakiinnitteisen maton kummaltakin puolelta.
- 2. Upota alusta veteen tai huuhtele se perusteellisesti.
- 3. Jätä matto tippumaan.
- 4. Levitä kostea matto puhtaalle alustalle ja asemoi se paikalleen.
- 5. Anna maton kuivua alustalla 24 tuntia ennen käyttöä.

#### Astian ominaisuudet

Suurin nopeus määräytyy astian tyypin mukaan, ja se määräytyy koon, muodon, pohjan muodon, materiaalin, pinnan epätasaisuuden ja puhtauden mukaan. Nämä parametrit tulee arvioida, jotta voidaan määrittää parhaiten käyttötarkoitukseen soveltuva matto ennen ravistimen käyttöä.

Käytä Taulukko 53 halutessasi määrittää, soveltuuko laboratoriolaitteistosi materiaali käytettäväksi liimakiinnitteisten mattojen kanssa.

Muovit/lasitavara	Tahmeusluokitus	Metalli	Tahmeusluokitus				
Lasi (borosilikaattilasi)	5	Ruostumaton teräs	5				
Polypropyleeni (PP)	3	Alumiini	5				
Polypropeenin kopolymeeri (PPCO)	3	Alumiini, eloksoitu	4				
Polyetyleeni (HPDE) (LPDE)	3						
Polykarbonaatti (PC)	4						
Polyeteenitereftalaatti (PET)	4						
Polyeteenitereftalaatin kopolyesteri (PETG)	4						
Polymetyylipenteeni (PMP)	3						
Polytetrafluorieteeni (PTFE)	0*						
Tetrafluorietyleeni- heksafluoripropeenin kopolymeeri (FEP)	0*						
Perfluorialkoksipolymeerit (PFA)	2						
Silikoni	1*						
Luokitukset: 0 = ei lainkaan tahmea; 5 = paras tahmeus / * Älä käytä astioita, joiden luokitus							

on 0 ja 1 (PTFE, FEP ja silikoni).

Taulukko 53: Astiamateriaalien yhteensopivuus

HUOMIO Suurin ravistelunopeus on saavutettavissa vain materiaaleilla, joiden luokitus on "5", ks. Taulukko 53.

HUOMIO Muut materiaalit kuin lasi tulee testata, ja sinun tulee itse määrittää, millaista materiaalia voidaan käyttää minkäkin kuormituksen, nopeuden, käyttöajan ja lämpötilan yhteydessä.
### Lasiastian (täyttöaste 30 %) nopeusalue ja käyttöaikaraja

Lasiastian nopeuden ja käyttöajan rajoitukset huoneenlämmössä (täyttöaste 30 %). Korkeampi lämpötila vähentää tahmeutta. Lämpötilan muuttaminen voi irrottaa astiat matosta. Aloita ravistelu akklimatisoiduilla astioilla.

	Mikrolevy / Deepwell-levy	Dekantterilasi						
	96 x 2 ml	25 ml	125 ml	250 ml	500 ml			
125 rpm	ei käyttöaikarajaa							
150 rpm	72 h	72 h	72 h	72 h	72 h			
200 rpm	72 h	72 h	72 h	72 h	72 h			
250 rpm	72 h							
300 rpm								
350 rpm								
400 rpm								

Taulukko 54: Kevyesti kiinnittyvä matto – nopeusalue ja käyttöaikaraja huoneenlämmössä

	Mikrolevy / Deepwell-levy	Erlenm	neyer						
	96 x 2 ml	25 ml	125 ml	250 ml	500 ml	1000 ml	2000 ml	3000 ml	5000 ml
125 rpm	ei käyttöaikarajaa	a							
150 rpm	72 h	72 h	72 h	72 h	72 h	72 h	72 h	72 h	72 h
200 rpm	72 h	72 h	72 h	72 h	72 h	72 h	72 h	72 h	72 h
250 rpm	72 h	72 h	72 h	72 h	72 h	72 h	72 h	48 h	24 h
300 rpm	72 h	72 h	72 h	72 h	72 h	72 h	48 h	24 h	
350 rpm	48 h	48 h	48 h	48 h	48 h	48 h	24 h		
400 rpm	24 h	24 h	24 h	24 h	24 h				

Taulukko 55: Tiukasti kiinnittyvä matto – nopeusalue ja käyttöaikaraja huoneenlämmössä

### <u>Valmistelu</u>

NOUDATA VAROVAISUUTTA Käytä aina suojalaseja ja -käsineitä käsitellessäsi lasitavaraa.

Ennen kuin laitat astiat liimakiinnitteiselle matolle:

- 1. Tarkista, ettei astioissa ole vaurioita. Älä käytä vahingoittuneita astioita.
- 2. Pyyhi astia puhtaalla ja kuivalla pyyhkeellä, jotta saat poistettua kaiken kosteuden.
- 3. Poista kaikki rasva ja muut epäpuhtaudet.
- 4. Tarkista, että astian pohja ja liimakiinnitteinen matto ovat kuivat ja puhtaat.

NOUDATA VAROVAISUUTTA Vaikka säiliön pohjassa olisi vain vähän näkyvää kalvoa tai yksittäinen vesipisara, säiliön kiinnitys voi löystyä ja säiliö voi irrota alustasta tietyn käyttöajan jälkeen.



Kuva 97: Liimakiinnitteinen matto: Pyyhi astia kuivaksi

5. Vähennä täyttömäärää tai sulje säiliöt, jotta nestettä eiroisku.

NOUDATA VAROVAISUUTTA Vaikka astia kiinnittyisikin kunnolla, yksikin tippa ulkopuolta pitkin valuvaa nestettä voi saada säiliön irtoamaan matosta tietyn käyttöajan jälkeen.

#### Astioiden asennus

- 1. Aseta astiat matolle ja paina niitä alaspäin.
- 2. Paina tai vedä varovasti, jotta jokainen säiliö tarttuu varmasti alustaan.

HUOMIO Pidä mielessäsi, että kondensoitunutta vettä voi muodostua alhaisessa lämpötilassa tai pitkäaikaisessa käytössä, mikä voi aiheuttaa astioiden irtoamisen vahingossa. Näytteen ja astian lämpötila ei saa poiketa ravistimen määritetystä lämpötila-alueesta. Lisätietoja ravistimen mahdollisesta lämpötila-alueesta on kohdassa "1. 1. Tekniset tiedot" sivulla 11.

### Astioiden irrotus

- 1. Paina näytön pysäytyspainiketta tai odota, kunnes käyttö on päättynyt.
- 2. Odota, että alusta pysähtyy kokonaan.
- Vedä astiaa sen kaulasta varovasti ja tasaisesti sivulle ja odota muutama sekunti. Jos astioiden pohjapinta-ala on suuri, niiden irtoaminen voi kestää muutaman sekunnin.



Kuva 98: Liimakiinnitteinen matto: Paina tai vedä astia varovasti irti

NOUDATA VAROVAISUUTTA Paina tai vedä varovasti äläkä käytä kovin paljon voimaa! Astiat saattavat irrota äkillisesti ja niiden sisältö voi läikkyä!

4. Jos astia ei irtoa, kaada vähän vettä sen alaosan ympärille.

Anna maton kuivua ennen kuin käytät sitä uudelleen.

#### Maton irrotus

1. Irrota liimakiinnitteinen matto nostamalla sitä toiselta puolelta ja vetämällä se hitaasti pois alustalta, ks. Kuva 99.



Kuva 99: Liimakiinnitteinen matto: Irrota matto nostamalla ja vetämällä se pois

#### Puhdistus ja desinfiointi

Jos tahmeus vähenee, puhdista matto.

NOUDATA VAROVAISUUTTA Älä käytä muita liuottimia, pesuaineita tai puhdistusaineita kuin mietoa astianpesuainetta, jossa ei ole ihonhoitoon tarkoitettuja lisäaineita. Suositeltava pitoisuus on enintään 2 ml pesuainetta 1 litrassa vettä.

#### Tarvikkeet

**HUOMIO** Maton suora kosketus pesuainetiivisteeseen vähentää tahmeutta. Tällöin pelkkä huuhtelu ei riitä. Jos käytät pesuainetta suoraan matolle, huuhtele matto, liota sitä vedessä 24 tuntia ja kuivaa sitä sitten 24 tunnin ajan, jotta liimautuvuus palaa täysin ennalleen.

HUOMIO Desinfioi ainoastaan 70-prosenttisella isopropanolilla tai 70-prosenttisella etanolilla.

#### Hieman värjäytynyt tai pölyyntynyt:

- 1. Pyyhi märällä tai desinfioi alustan matot päältä.
- 2. Kostuta matto.
- 3. Pyyhi jäämät pois sienellä.
- 4. Pyyhi puhtaaseen veteen kostutetulla sienellä.
- 5. Desinfioi tarvittaessa.
- 6. Anna kuivua 24 tuntia.

#### Erittäin likainen:

- 1. Irrota matto alustasta, jotta voit puhdistaa maton kokonaan.
- 2. Liota mattoa veteen laimennetussa astianpesusaippuassa.
- 3. Poista kaikki jäämät hankaamalla alustaa huolellisesti sienellä.
- 4. Huuhtele alusta puhdistuksen jälkeen riittävällä vedellä.
- 5. Desinfioi tarvittaessa.
- 6. Anna kuivua 24 tuntia.
- 7. Suorita asennus, ks. kohtaa "Kuivamaton asennus" sivulla 107 tai "Märkämaton asennus" sivulla 107.

### 3. 3. 12. Kaasunjakoputkiston asennus

Kaasunjakoputkisto mahdollistaa inertin kaasun, kuten typen tai hiilidioksidin, injektoinnin useisiin astioihin ravistimen käytön aikana (katso luvusta "1. 2. 12. Yleiset lisätarvikkeet" sivulla 42 tarvittavat tilaustiedot).

Kaasunjakoputkisto on asennettu tehtaalla Solaris 2000 I/R - ja 4000 I/R -ravistimiin, ja ne ovat saatavilla myös asiakkaan asennettavana lisätarvikkeena, jos niitä ei ole alun perin asennettu yksikköön. Kaikissa Solaris 2000 I/R - ja 4000 I/R -ravistimissa on käyttöportti kummassakin sivupaneelissa (katso luvusta "1. 4. Tuotekatsaus" sivulla 44 tarvittavat sijaintitiedot) laboratoriokaasun syöttöletkun kautta imusarjaan tapahtuvaa syöttöä varten. Kaasun jakoputkistossa on astioille yksi kaasun syöttöliitin ja kahdeksan poistoliitintä.



Kuva 100: Solaris 4000 R -ravistimeen asennettu kaasunjakoputkisto

#### Suorita asennus seuraavasti:

- 1. Kiinnitä kaasunjakoputkisto ② tukevasti ravistimen kammion takaseinään toimitukseen sisältyvillä ruuveilla ⑤.
- Irrota muovitulppa käyttöportin läpivientikumin ulkopuolelta (katso luvusta "1. 4. Tuotekatsaus" sivulla 44 tiedot sijainnista).
- 3. Tee veitsellä ristikkäinen viilto ravistimen kammion sisäpuolella olevan läpivientikumin kalvoon.
- 4. Leikkaa pala joustavaa putkea sopivan pituiseksi laboratoriokaasun syöttöputkeksi ①.
- 5. Kiinnitä laboratoriokaasun syöttöputki jakoputkiston ① kynnelliseen syöttöliittimeen ④.

- 6. Varmista, että laboratoriokaasun syöttöletku ① pysyy irti kannen vivuista, ja työnnä se käyttöaukon kumisen läpivientikukin läpi ulos.
- 7. Kiinnitä laboratoriokaasun syöttöletku ① kaasulähteeseen.
- Leikkaa 8 putkenpalaa ja kiinnitä ne jakoputkiston ulostuloihin ③. Jos astioita on yli 8, käytä Y-liittimiä. Jos astioita on vähemmän kuin 8, purista käyttämättömät letkut kiinni kaasuhävikin vähentämiseksi.
- 9. Työnnä putket astioihin sopiva tulppia ja tiivisteitä käyttämällä.

### 3.4. Lastaus ja normaali käyttö



# Tulipalovaara johtuen käynnistyneistä kemiallisista reaktioista.

VAROITUS

Älä käytä ravistinta nopeuksilla, jotka heittävät ulos astioiden sisällön.

Lisää nopeutta hitaasti. Kokeile vedellä ennen kemikaaleja.



# Turvallisuus voi vaarantua väärästä lastaamisesta ja rikkonaisista apuvälineistä.



- Varmista, että lasti (apuvälineet ja näytteet) on sijoitettu symmetrisesti alustan keskukseen nähden. Kun käytät dual stack -alustaa, varmista, että alemmalla alustalla on suurempi kuorma kun ylemmällä.
- Älä ylikuormita ravistinta. Katso enimmäiskuormitus kohdasta "Tekniset tiedot" sivulla 11. Kuorma sisältää alustan painon, ravistimen päälle asetettujen apuvälineiden ja näytteiden painon.
- Huolehdi, että apuvälineet on asennettu oikein ennen kuin käytät ravistinta. Noudata ohjeita osiossa "Tarvikkeet" sivulla 88.



#### Sulkeutuva kansi voi aiheuttaa sormi- tai käsivamman.

 Ravistimen kansi voi pudota vahingossa ja kiilata tai puristaa sormia tai käsiä ja aiheuttaa näin vammoja. Näin voi käydä, kun kannen jousi on viallinen eikä pysty pitämään kantta täysin auki-asennossa. Älä käytä ravistinta, jonka jousi on viallinen. Ota yhteys Thermo Fisher Scientificin asiakaspalveluun ja vaihdata kannen jousi välittömästi.



### Lataa

Huolehdi, että kuorma, mukaan lukien apuvälineet ja näytteet, on asetettu symmetrisesti alustan keskelle. Kun käytät dual stack-alustaa etkä käytä koko kapasiteettia: Jotta saisit parhaan tuloksen, lastaa mieluiten ala-alusta ja lastaa symmetrisesti alustan keskelle.

Varmista lämpötilaohjattua ravistinta käyttäessäsi, että kuorma on asennossa, joka jättää riittävästi tilaa kannen sulkemiseen. Varmista myös, että tarvittava liike on mahdollinen kansi suljettuna, jotta törmäyksiä tai vaurioita ei synny.



Kuva 101: Esimerkki hyvin lastatusta alustasta

												ā	Ā
(•	٠	٠	٠		*		*	*	÷	٠	-	•	$\sim$
•	٠	٠	*	٠	*	٠	*	*	٠	٠	. "	÷	·(·)) · ·
•	٠	٠	٠	*	*	٠	*	*	٠	*	*	۴	
•			*	٠	*		*			*	*	*	· 🉈 🕄 🕩
•	٠	*	٠	*	٠	*	*	*		*	٠	٠	<b>((·)</b> ) <sup>1</sup> <sup>∞</sup>
•			*		*			·	·	٠		٠	
•	•	,	•	*	•	*	*	*		*	٠	٠	· 🙈 📢 🕩
•		*	*			*	*	*		*	*	*	
•	٠	٠	•	•	٠	•	٠		٠	٠	۰	٠	
•				*		*		*		٠	*	e	· 🙈 🌔 🕽
Ŀ	•	•	•	•	•	•	*	٠	٠	٠	٠	٠	·(·)• · · ·

Kuva 102: Esimerkki huonosti lastatusta alustasta

HUOMIO Maksimilasti sisältää alustan painon, apuvälineet ja näytteet, jotka on asennettu ravistimeen. Katso kohdasta "1. 1. Tekniset tiedot" sivulla 11 tiedot maksimikuormituksesta.

### Normaali käyttö

Tiedot nopeuskuorman rajoituksista annetaan vain ohjeellisina tuotteen käyttöä varten. Tasainen käynti matalalla tärinätasolla tai kokonaan ilman tärinää riippuu useista tekijöistä kuten täyttöasteesta, pidikkeiden tyypistä ja kunnosta, astian tyypistä, astioiden järjestelystä alustalla, pinnasta, jolle ravistin on asetettu, ja valitusta nopeudesta.

Kun asetat uuden tai tuntemattoman yhdistelmän tai kun joku näistä parametreista muuttuu: Lisää nopeutta yli 175 rpm hitaasti ja seuraa ei-haluttua lastin tärinää tai ravistimen siirtymistä. On sinun vastuullasi käyttää ravistinta turvallisesti.

Solaris-ravistimet on suunniteltu erityisesti vähän lämpöä ja vähän tärinää tuottaviksi, joten ne sopivat erinomaisesti käytettäväksi useissa laboratoriolaitteissa, kuten ilmastoiduissa kammioissa, inkubaattoreissa ja laboratoriojääkaapeissa. Erityisten ympäristöolosuhteiden, kuten hyllyjen ja kannattimien vakauden, vaihtelun vuoksi käyttäjä on vastuussa turvallisesta käytöstä, kun laitetta käytetään jossakin laboratoriolaitteessa. Lisää nopeutta hitaasti ja tarkista, ettei laite liiku tahattomasti.

#### Turvallinen nopeus

Solaris-ravistimien kuormituksesta riippumaton turvanopeus näkyy Taulukko 56. Kuorma sisältää alustan painon, ravistimen päälle asetettujen apuvälineiden ja näytteiden painon. Katso kohdasta "Tekniset tiedot" sivulla 11 tiedot maksimikuormituksesta.

Malli	Turvallinen enimmäisnopeus
Solaris 2000	175 rpm
Solaris 4000	175 rpm
Solaris 2000 I	250 rpm
Solaris 2000 R	250 rpm
Solaris 4000 I	250 rpm
Solaris 4000 R	250 rpm

Taulukko 56: Turvallinen enimmäisnopeus

#### Korkein nopeus

Korkeimmalla 525 rpm nopeudella Solaris-ravistimia voidaan käyttää rajallisella kuormituksella, ks. Taulukko 57. Kuorma sisältää alustan painon, ravistimen päälle asetettujen apuvälineiden ja näytteiden painon. Huomioi, että käytettyjen puristimien ja astioiden nopeuskapasiteetti ei välttämättä sovellu kyseiselle nopeudelle.

Malli	Turvalliset enimmäisnopeudet
Solaris 2000	3 kg
Solaris 4000	10 kg
Solaris 2000 I	4 kg
Solaris 2000 R	4 kg
Solaris 4000 I	8 kg
Solaris 4000 R	8 kg

Taulukko 57: Enimmäiskuormitus suurimmalla nopeudella

HUOMIO Aina kun nopeus lisääntyy, puristimet alkavat aueta ja lisätä ylimääräistä tilapäistä tärinää järjestelmään johtuen astioiden suuremmasta liikkeestä.

#### Paino ja nopeus -käyrät

Seuraavien paino ja nopeus -taulukkojen avulla voit määrittää parhaan nopeuskuormitus -yhdistelmän normaalikäyttöösi. Vihreä alue näyttää nopeuslastiyhdistelmät, jotka eivät aiheuta tärinää tai aiheuttavat vain pientä tärinää sekoituksen aikana. Punainen alue näyttää yhdistelmät, jotka voivat aiheuttaa voimakasta tärinää ravistelun aikana ja voivat johtaa ravistimen tafattomaan liikkumiseen. Koska nämä kaaviot ovat vain suuntaa-antavia, sinun on oltava tarkkana, kun nopeus-lastiyhdistelmä lähenee punaista aluetta.

Huomioi, että käytöllesi ominaiset olosuhtet voivat aiheuttaa ei-haluttua lasti- tai instrumenttikäytöstä, ennen kuin ne saavuttavat näytetyn rajan. Lisää nopeutta vaiheittain nähdäksesi määrätyn lastin käyttäytymisen. Kuorma sisältää alustan painon, ravistimen päälle asetettujen apuvälineiden ja näytteiden painon.

HUOMIO Seuraavien kaavioiden sininen vaakaviiva osoittaa kunkin alustan tyhjäpainon.

#### Solaris 2000



Kuva 103: Solaris 2000 – Normaali käyttö



#### Solaris 4000



Lastaus ja normaali käyttö



Kuva 105: Solaris 2000 I / 2000 R - Normaali käyttö



#### Solaris 4000 I / 4000 R

Kuva 106: Solaris 4000 I / 4000 R - Normaali käyttö

# 4. Huolto ja kunnossapito



VAROITUS

#### Vaarallisten aineiden käsittelyn aiheuttama riski

- Mmitä tahansa vaarallisia materiaaleja sekoitettaessa on huomioitava Maailman terveysjärjestön (WHO) "Laboratorion bioturvallisuuskäsikirja" ja paikalliset säädökset. Kun sekoitetaan mikrobiologisia näytteitä riskiryhmässä II (Maailman terveysjärjestön (WHO) "Laboratorion bioturvallisuuskäsikirjan" mukaan), on käytettävä aerosolivarmoja biologisia tiivisteitä. Katso internetistä sivua World Health Organization (www. who.int) löytääksesi "Laboratory Biosafety Manual'in". Korkeamman riskin ryhmien materiaaleille on käytettävä erikoisturvatoimenpiteitä.
- Jos toksiinit tai patogeeniset aineet ovat kontaminoineet ravistimen tai sen osia, tarvitaan desinfiointitoimia ("Dekontaminaatio" sivulla 124; "Desinfiointi" sivulla 124).
- Jos vaarallinen tilanne sattuu, käännä virta pois ravistimesta ja poistu välittömästi alueelta.



VAROITUS

#### Vauriot terveydelle tartuntavaarallisista aineista

Jos nesteitä tai muita aineita vahingossa läikkyy alustan alle, sammuta ravistin välittömästi, irrota se pistorasiasta ja poista alusta ("Alustat" sivulla 18).

Puhdista läikkynyt materiaali seuraten tavanomaisia laboratoriotoimenpiteitä. Käytä kunnollisia henkilökohtaisia suojaimia.

Kaikki sisäiset säädöt ja korjaukset on tehtävä valtuutetun huoltoteknikon toimesta. Käyttäjä ei saa avata ravistimen koteloa.

Noudata kaikkia kyseisten apuvälineiden mukana toimitettuja tuotetietoja, joissa kerrotaan tarkasti, kuinka tuote huolletaan ja puhdistetaan oikein. Käytä tämän kappaleen sisältämiä tietoja vain suuntaa antavina ohjeina.

### 4.1. Perusteita

Henkilökohtaisen, ympäristöllisen ja materiaaliturvallisuuden takia sinun tulee puhdistaa, ja jos tarpeellista, desinfioida ravistin ja apuvälineet säännöllisesti.

Thermo Fisher Scientific suosittelee laboratorioravistimen puhdistusta ja manuaalista desinfiointia vähintään kerran kuussa. Normaali huoneilma sisältää tuhansia kierteleviä mikro-organismeja, jotka voivat pesiytyä ravistimeen vaarantaen näin viljelmäsi.

Thermo Fisher Scientific suosittelee käyttämään 70 % etanolia, tai 70 % isopropanolia tai korkeintaan 10 % ammoniumpohjaista desinfiointiainetta.

HUOMIO Mekanismi voi vaurioitua nesteiden sisään menosta. Älä päästä nesteitä, erityisesti orgaanisia liuottimia, kontaktiin mekanismin tai sen laakeroinnin kanssa. Orgaaniset liuottimet hajottavat rasvan mekanismin laakeroinnissa. Kun nesteitä tihkuu ravistelijan sellaisiin osiin, joita ei voi puhdistaa, älä käytä ravistinta ja ota yhteys huoltoon.

HUOMIO Ei-hyväksytyt toimenpiteet tai aineet voivat heikentää ravistimen materiaaleja ja johtaa käyttöhäiriöön. Älä käytä mitään puhdistus- tai dekontaminaatiotoimenpidettä, jollet ole täysin varma aiotun toimenpiteen turvallisuudesta laitteistolle. Käytä vain puhdistusaineita, jotka eivät vahingoita laitteistoa. Jos epäilet yhteensopivuutta ota yhteys puhdistusaineiden valmistajaan.

**HUOMIO** Huolehdi lämpötilaohjattuja ravistimia käyttäessäsi siitä, että jätät kannen auki, jotta puhdistusaineet ja vesi pääsevät haihtumaan.

- Irrota verkkojohto ennen puhdistusta, desinfiointiaa tai dekontaminaatiota.
- Poista asennetut apuvälineet ja alusta(t) ravistimesta ennen puhdistusta, desinfiointia ja dekontaminaatiota.
- Käytä lämmintä vettä miedolla puhdistusaineella ja pehmeää liinaa materiaalien puhdistuksen. Ongelmatapauksissa ota yhteys Thermo Fisher Scientificiin. Huuhtele puhtaalla vedellä ja kuivaa läpikotaisin.
- Älä koskaan käytä syövyttäviä puhdistusaineita, kuten fosforihappoa, valkaisuaineita tai hankausjauhetta.
- Käytä vain desinfektioaineita, joiden pH on 6–8.
- Puhdista läikkeet välittömästi nukattomalla liinalla, joka on kostutettu ei-syövyttävällä puhdistusaineella, puhdistusaineen valmistajan ohjeiden mukaan.
- Läikkeet voivat valua alustan alle. Jos alustan alle läikkyy jotakin, irrota alusta ja poista kaikki läikkynyt aine alustan alla olevalta keräystasolta. Desinfioi tai dekontaminoi tarvittaessa.
- Tarkista varistimen osat ja poista kaikki läikkynyt kasvualusta tai jäänteet.

### Apuvälineiden tarkistus

HUOMIO Älä käytä mitään ravistinta tai apuvälineitä, joissa on vaurion merkkejä. Suosittelemme, että tutkitutat apuvälineet säännöllisesti osana rutiinihuoltoa varmistaaksesi turvallisuuden.

Kun apuvälineet on läpikotaisin puhdistettu, täytyy ne tutkia vaurioiden, kulumisen ja korroosion varalta.

#### <u>Metalliosat</u>

Jos vaurioita löytyy, kuten korroosiota, kulumista tai lohkeamia, apuvälineet täytyy poistaa välittömästi käytöstä.

#### <u>Muoviosat</u>

Tarkistus säröjen, haalistumisen, pintavikojen tai lohkeamien varalta. Vian löytyessä tutkittava osa on poistettava välittömästi käytöstä.

#### Liimakiinnitteiset matot

Tarkista ennen jokaista käyttökertaa, että liimakiinnitteisissä matoissa ei ole kulumisen tai vaurioitumisen merkkejä, kuten säröilyä, värinmuutoksia, kellastumista, haurastumista, muodonmuutoksia, pinnan hankaumia tai kemiallisia vaurioita. Hävitä välittömästi liimakiinnitteiset matot, joissa on merkkejä kulumisesta tai vaurioista.

### 4.2. Puhdistus

HUOMIO Ennen kuin käytät mitään puhdistusmetodeja, käyttäjän tulee tarkistaa valmistajalta, etteivät puhdistusaineet tai aiottu metodi vahingoita laitetta.

HUOMIO Mekanismi voi vaurioitua nesteiden sisään menosta. Älä päästä nesteitä, erityisesti orgaanisia liuotteita, mekanismiin tai mekanismin laakerointiin. Orgaaniset liuottimet hajottavat rasvan mekanismin laakeroinnissa. Pese yksikön ulkopuoli pehmeällä liinalla käyttäen mietoa saippuavesiliuosta, huuhdo puhtaalla vedellä ja kuivaa perusteellisesti.

Katso kohdasta "Perusteita" sivulla 122 tiedot ravistimen ja käytettyjen tarvikkeiden oikeasta puhdistuksesta.

**HUOMIO** Huolehdi lämpötilaohjattuja ravistimia käyttäessäsi siitä, että jätät kannen auki, jotta puhdistusaineet ja vesi pääsevät haihtumaan.

### Kosketusnäyttö

Puhdistaaksesi kosketusnäytön:

- 1. Irrota sähkökosketin seinästä.
- 2. Puhdista näyttö mikrokuituliinaa käyttäen.
- 3. Jos tarpeen, kostuta mikrokuituliina vedellä ja pyyhi kosketusnäyttö uudelleen.
- 4. Anna kuivua.

### 4.3. Desinfiointi

VAROITUS Riski vaarallisten aineiden käsittelystä. Älä koske infektoituneisiin osiin. Vaarallinen infektio on mahdollinen kun kosketaan kontaminoituneisiin osiin. Tartuntavaarallinen materiaali voi päästä ravistimeen, kun astia rikkoontuu, tai läikkymisen seurauksena. Kontaminaation sattuessa varmistu, ettei kukaan joudu vaaraan. Desinfioi kyseiset osat välittömästi.

HUOMIO Tarvikkeet voivat vaurioitua vääristä desinfiointimetodeista tai-aineista. Varmistu, ettei desinfektioaine tai-metodi vahingoita laitteita. Jos olet epävarma, ota yhteyttä desinfektioaineen valmistajaan. Seuraa desinfektioaineiden turvatoimenpiteitä ja käsittelyohjeita.

- 1. Pyyhi kaikki osat ja alueet 70 % etanolilla halutun desinfiointiasteen mukaan. Älä kastele mitään alueita, joissa on paljaana elektroniikkaa.
- 2. Anna kuivua.

HUOMIO Huolehdi lämpötilaohjattuja ravistimia käyttäessäsi siitä, että jätät kannen auki, jotta puhdistusaineet ja vesi pääsevät haihtumaan.

Olet vastuussa siitä, että haluamasi desinfiointiaste saavutetaan asianmukaisesti.

### 4.4. Dekontaminaatio

VAROITUS Riski vaarallisten aineiden käsittelystä. Älä koske kontaminoituneisiin osiin. Altistus kontaminaatiolle on mahdollinen kun, kontaminoituneisiin osiin kosketaan. Kontaminoitunut materiaali voi päästä ravistimeen, kun putki menee rikki, tai läikkymisen seurauksena. Kontaminaation sattuessa varmistu, ettei kukaan joudu vaaraan. Dekontaminoi kyseiset osat välittömästi.

HUOMIO Laitteisto voi vaurioitua sopimattomista dekontaminaatiometodeista tai-aineista. Varmistu, ettei dekontaminaatioaine tai -menetelmä vahingoita laitteita. Epäselvissä tapauksissa ota yhteys dekontaminaatioaineen valmistajaan. Seuraa desinfektioaineiden turvatoimenpiteitä ja käsittelyohjeita

Thermo Fisher Scientific suosittelee seuraavaa metodia.

- 1. Pyyhi kaikki osat ja alueet 70 % etanolilla. Älä kastele mitään alueita, joissa on paljaana elektroniikkaa.
- 2. Anna kuivua.

HUOMIO Huolehdi lämpötilaohjattuja ravistimia käyttäessäsi siitä, että jätät kannen auki, jotta puhdistusaineet ja vesi pääsevät haihtumaan.

Olet vastuussa siitä, että haluttu dekontaminaatioaste saavutetaan.

### 4.5. Autoklavointi

HUOMIO Älä koskaan ylitä autoklavoitaessa sallittua lämpötilaa tai aikaa. Höyryssä ei saa olla mitään kemikaaleja/lisäaineita.

Ennen autoklavointia:

- Pura apuvälineet.
- Huuhdo kunnolla pois kaikki kemikaali- ja puhdistusainejäämät. Tee lopullinen huuhtelu tislatulla vedellä.

Tästä taulukosta näet autoklavointimahdollisuudet:

Esine	Autokla- voitava	Tarkemmat tiedot
Ravistin	Ei	-
Alusta	Kyllä	121 °C, 20 min
Puristin (ilman jousiputkia)	Kyllä	121 °C, 20 min
Ruuvitaltta	Ei	-
Mikrolevyn / Deepwell-levyn pidike	Kyllä	121 °C, 20 min
Kiinnittyvä matto	Ei	-
Dekantterilasiteline	Kyllä	121 °C, 20 min
Erotussuppilon pidike	Kyllä	121 °C, 20 min
Koeputkiteline – metalliosat	Kyllä	121 °C, 20 min
Koeputkiteline – muoviosat	Ei	-
Koeputkiteline – vaahtomuovityynyt	Ei	-
Säädettävä astianpidike – metalliosat	Kyllä	121 °C, 20 min
Säädettävä astianpidike – liukumaton matto	Ei	-
Tarviketaso – metallialusta	Kyllä	121 °C, 20 min
Tarviketaso – kumimatto	Ei	-
Kallistuskulmaltaan säädettävän koeputkitelineen pidike – metalliosat	Kyllä	121 °C, 20 min
Kallistuskulmaltaan säädettävän koeputkitelineen pidike – muoviosat	Ei	-
Kallistuskulmaltaan säädettävän koeputkitelineen pidike – vaahtomuovityynyt	Ei	-

Taulukko 58: Materiaalien autoklavoitavuus

Pidä huolta, että vaatimusten mukainen steriliteetti saavutetaan.

Ulkonäkö ja väri voivat muuttua hiukan autoklavoitaessa.

### 4.6. Lämpötilan kalibrointi

Lämpötilaohjatun ravistimen lämpötilan asetusarvo voidaan kalibroida tunnetulla tarkkuudella mittaavan lämpötilan mittauslaitteen avulla, jotta voidaan varmistaa inkuboidun tai jäähdytetyn ravisteluprosessin täydellinen toistettavuus.

### 4. 6. 1. Lämpötilan kalibroinnin parhaat käytännöt

Noudata ohjeita tarkasti ennen ravistimen lämpötilan kalibrointia ja sen aikana. Mahdolliset häiriöt tai poikkeamat suositelluista parhaista käytännöistä voivat estää näytteen lämpötilan stabiloitumisen ja aiheuttaa kalibroinnin epäonnistumisen.

Parhaita käytäntöjä ovat mm. seuraavat:

- 250 ml:n mittapullon käyttö.
- Pullon täyttö täsmälleen 100 ml:lla vettä.
- Sulje kansi kokonaan ja pidä se suljettuna koko kalibroinnin ajan.
- Anna näytteen lämpötilan stabiloitua riittävän kauan, jotta lämpötila stabiloituu täysin.
- Älä anna ympäristön olosuhteiden, kuten lämpötilan ja kosteuden, muuttua kalibroinnin aikana. Pienetkin muutokset, kuten oven tai ikkunan avaaminen tai ilmastointijärjestelmän asetusten muuttaminen, voivat vaarantaa käynnissä olevan kalibrointiprosessin tuloksen.
- Ravistin kannattaa kalibroida vuosittain.

### 4. 6. 2. Lämpötilan kalibrointimenettely

Suorita kalibrointi seuraavasti:

- 1. Voit käynnistää kalibroinnin kolmella eri tavalla:
  - » Kun alkukäynnistys on suoritettu kohdassa "Ensimmäinen käynnistys" sivulla 54 kuvatulla tavalla, voit suorittaa lämpötilan kalibroinnin sen sijaan, että poistuisit alkukäynnistyksestä.
  - » Ravistin näyttää kalibrointimuistutuksen vuosittain.
  - » Tämän jälkeen voit käynnistää kalibroinnin napauttamalla valikkorivin Asetukset -kuvaketta ja painamalla sen jälkeen Kalibrointi -painiketta alempana Lämpötilan kalibrointi -näytöllä.
  - » Napauttamalla Asetukset-kuvaketta valikkorivillä ja sen jälkeen Kalibrointi-painiketta, voit käynnistää prosessin milloin tahansa tarvittaessa.

HUOMIO Kalibrointimuistutus voidaan poistaa käytöstä **Poista kalibrointi-ilmoitukset** käytöstä -valintaruudulla Asetukset -näytöstä.

2. Odota, että Lämpötilan kalibrointi -näyttö, ks. Kuva 107, tulee näkyviin.



Kuva 107: Kalibrointi: Aloitusnäyttö

3. Napauttamalla Aloita kalibrointi -painiketta pääset aloittamaan kalibroinnin.



Kuva 108: Kalibrointi: Erlenmeyer-pullon täyttökehote

- 4. Noudata näyttöön tulevia ohjeita: Täytä 250 ml:n erlenmeyerpullo100 ml:lla vettä.
- 5. Aseta pullo pidikkeeseen ja kiinnitä pidike suunnilleen ravistelualustan keskelle.
- Ripusta mittapulloon tunnetulla tarkkuudella mittaavan lämpötilan mittauslaitteen anturi. Anturi tulee upottaa nesteeseen, mutta se ei saa koskettaa pullon pohjaa tai sivuja.
- 7. Sulje pullo, jotta nestettä ei läiky.
- 8. Sulje ravistelualustan päällä oleva kansi.

9. Napauttamalla **Poista**-painiketta pääset jatkamaan.



Kuva 109: Kalibrointi: Aloita kalibrointi -kehote

HUOMIO Ravistin aloittaa ravistelun heti, kun napautat Käynnistä-painiketta.

 Noudata näyttöön tulevia ohjeita: Käytä esiasetettua lämpötilaa napauttamalla Käynnistä painiketta tai napauta Lämpötilan asetuspiste -kenttää ja syötä haluamasi lämpötila ennen kuin napautat Käynnistä-painiketta.

Näytössä näkyy, miten ravistimen kammio lämpiää (tai jäähtyy) valittuun kalibrointilämpötilaan.



Kuva 110: Kalibrointi: Kammion stabiloituminen kalibrointilämpötilaan

11. Odota vähintään kaksi ja puoli tuntia (150 minuuttia), jotta kammion lämpötilan mittaus saavuttaa vakaan tilan.

Vakaan tilan ilmaisee ilmoitus **Kammion lämpötila on stabiloitunut**, ks. Kuva 111 alla.



Kuva 111: Kalibrointi: Kammion stabiloituminen kalibrointilämpötilaan onnistui

- 12. Noudata näyttöön tulevia ohjeita:
  - » Seuraa termoelementin lämpötilalukemaa säännöllisesti sen jälkeen, kun Kuva 111 näyttö on tullut näkyviin. Jos katsot lämpötilan olevan vakaa, voit joko napauttaa **Muokkaa**-painiketta ja korjata sitä tai **Valmis**painiketta napauttamalla viimeistellä kalibroinnin.



Kuva 112: Kalibrointi: Ravistimen näyttämä oikea lämpötila

- 13. Korjaa ravistimen näytössä näkyvä lämpötila valitsinkiekolla tarkan mittauslaitteen lukeman mukaiseksi.
- 14. Napauttamalla Tallenna painiketta saat muutoksesi tallennetuiksi.

15. Lämpötilan kalibrointi valmis -viesti ilmaisee, että prosessin suoritus onnistui.

<	Shake Lab 1 11:22:24AM Temp Calibration	
¢. (	Step 5 of 5	
邕	Temperature calibration complete.	
Ē		
â.		
	Re-Calibrate	<b>(</b> )

Kuva 113: Kalibrointi: Lämpötilan kalibrointi onnistui

Napauttamalla Kalibroi uudelleen voit toistaa kalibrointiprosessin.

### 4. 6. 3. Lämpötilan kalibroinnin tarkistus

Voit tarkistaa nykyisen kalibrointilämpötilan milloin tahansa:

- 1. Napauttamalla **Settings**-kuvaketta valikkopalkissa ja sn jälkeen **Kalibrointi**painiketta alempana näkyviin tulevalla **Lämpötilan kalibrointi** -näytöllä
- 2. Lämpötilan kalibrointi -näyttö, ks. Kuva 107, tulee näkyviin. Tässä näytössä näkyy senhetkinen Ravistimen kalibroitu lämpötila.

HUOMIO Lämpötilaohjatuissa ravistimissa voit myös etsiä kalibrointilämpötilaa näytössä, jossa lämpötila on määritetty. Esimerkki tästä on osiossa "Lämpötilan määritys" sivulla 64.

### 4.7. Laiteohjelmiston asennus

Kun napautat Laiteohjelmisto-painiketta näytössä Tiedostot ja tiedot (ks. "Tiedostot ja info" sivulla 77), sinua pyydetään antamaan järjestelmävalvojan salasana, jonka jälkeen näkyviin tulee Laiteohjelmisto-näyttö, ks. Kuva 114. Tässä näytössä voit asentaa uuden laiteohjelmiston USB-muistitikulta.

Voit saada huoltosopimuksen mukaisen uuden laiteohjelmiston latauslinkin huoltoteknikolta. Asenna laiteohjelmisto vain valtuutetuista lähteistä. Muista kuin valtuutetuista lähteistä peräisin olevat ohjelmistot voivat mitätöidä Thermofisher Scientificin myöntämän takuun.

- 1. Lataa ZIP-tiedosto huoltoteknikon sinulle toimittamasta linkistä.
- 2. Pura ZIP-tiedosto tyhjään USB-asemaan.

HUOMIO Varmista, että USB-asemassa on riittävästi tilaa ZIP-tiedoston purkamiseen.

- Jos laite on edelleen käynnissä, pysäytä ravistin painamalla Pysäytys-painiketta näyttöikkunassa.
- 4. Kuittaa mahdolliset hälytykset, ks. osiota "Tila" sivulla 66.

HUOMIO Jos yrität suorittaa päivityksen ravistimen ollessa käynnissä, näyttöön tulee ilmoitus, että **Ravistimen päivitys ei ole sallittu ravistimen ollessa käynnissä.** Ilmoituksessa on **Pysäytä ravistin** -painike, jolla voit poistua näytöstä ja suorittaa päivityksen.

- 5. Napauta Tiedostot ja tiedot-kuvaketta navigointipalkissa.
- 6. Vieritä **Tiedostot ja tiedot** -näyttö alalaitaansa, jotta saat näkyviin **Laiteohjelmisto**-painikkeen.
- 7. Napauttamalla painimetta **Laiteohjelmisto** voit avata Laiteohjelmisto-näytön, jossa näkyvät tähän mennessä asennetut laiteohjelmistoversiot.
- 8. Liitä USB-asema johonkin ravistimen USB-porttiin. USB-porttien sijainti on kuvattu osiossa "Tuotekatsaus" sivulla 44.
- 9. Anna järjestelmänvalvojan salasana.



Kuva 114: Laiteohjelmisto-näyttö

10. Napauta Päivitä laiteohjelmisto -painiketta.

HUOMIO Jos ravistin ei löydä USB-asemaa, sinua kehotetaan **asettamaan sopivaan porttiin USB-muistitikku, jossa on laiteohjelmisto, jonka haluat asentaa päivityksen aloittamiseksi.** Aseta USB-muisti paikalleen nyt tai tarkista, että se on kunnolla kiinni USB-portissa ja korjaa näin ongelma.  Jos laiteohjelmiston päivitys on tarjolla, Kuva 115 eritelty luettelo tulee näkyviin Laiteohjelmiston päivitys -näyttöön. Napauta laiteohjelmistokohdetta, jonka haluat päivittää.

Firmware Upgrade	×	Firmware Upgrade	×
Select firmware to upgrade: H.M.I. Main Controller Parameters		Select firmware to upgrade: H.H.L Main Controller Parameters	
Select		Select	

Kuva 115: Valitse asennettava laiteohjelmisto

HUOMIO Ajan tasalla olevat laiteohjelmistokohteet näkyvät harmaina eikä niitä voi valita. Jos päivitettäviä kohteita ei ole saatavilla, näyttöön tulee ilmoitus, että **Ohjelmistoon ei tällä** hetkellä ole saatavana uusia päivityksiä.

- 12. Napauta Valitse.
- Jos muistitikussa on useampi kuin yksi laiteohjelmistoversio, sinua pyydetään valitsemaan jokin niistä. Valitse laiteohjelmistoversio napauttamalla ja napauta sn jälkeen Valitse.



Kuva 116: Valitse asennettava laiteohjelmistoversio

HUOMIO Jos et ole varma asennettavasta versiosta, ota yhteyttä latauslinkin toimittaneeseen huoltoteknikkoon.

14. Näkyviin tulee luettelo päivitettävistä kohteista. Napauttamalla **OK** voit aloittaa päivityksen.



Kuva 117: Päivitettävien kohteiden luettelo

**HUOMIO** Graafinen käyttöliittymä hoitaa keskinäiset riippuvuudet puolestasi: Valitse yksilaiteohjelmistokohde päivitettäväksi, ja ohjelmisto lisää kaikki tarvittavat kohteet automaattisesti asennukseen.

HUOMIO Jos valitset Kuva 115 luettelosta H.M.I. -vaihtoehdon, koska haluat päivittää ainoastaan graafisen käyttöliittymän laiteohjelman, graafisen käyttöliittymän näyttö ei vastaa kosketusnäytön aktiivisen alueen muihin napautuksiin jopa minuuttiin laiteohjelmistopäivityksen aloittamisen jälkeen. <u>Älä</u> sammuta ravistinta tänä aikana, muussa tapauksessa päivitys ei onnistu. Ravistin käynnistyy uudelleen lyhyen ajan kuluttua ja on taas valmis käytettäväksi.

**HUOMIO** Jos valitset Kuva 115 luettelosta päivitettäväksi <u>useita</u> laiteohjelmistokomponentteja kerralla, ravistinta <u>ei saa</u> sammuttaa missään vaiheessa päivitysprosessin aikana.

15. Näyttöön tulee erilaisia viestejä, jotka kertovat asennuksen edistymisestä, ks. Kuva 118.



Kuva 118: Laiteohjelmistopäivityksen asennusaikaiset ilmoitukset

16. Kun sinua kehotetaan käynnistämään ravistin uudelleen, katkaise laitteesta virta ja käännä se uudelleen päälle virtakytkimellä.

### 4.8. Alustan vaihto

Alustan asennuksen perusmenettely on kuvattu luvussa "3. 3. 1. Alustan asennus" sivulla 89. Täydellinen luettelo kunkin ravistimen varaosa-alustoista on osiossa "1. 2. 1. Alustat" sivulla 18.

HUOMIO Lämpötilaohjattuja ravistinmalleja voidaan käyttää vain tehdasasenteisten alustojen kanssa.

### 4.9. Kunnossapito

Thermo Fisher Scientific suosittelee ravistimen ja apuvälineiden huoltoa vuosittain valtuutetun teknikon toimesta. Huoltoteknikko tarkistaa seuraavat:

- sähkölaitteet
- käyttöpaikan sopivuus
- turvajärjestelmä
- käytetyt apuvälineet
- puristimien ja alustojen sekä muiden apuvälineiden kiinnitys ravistimeen

Ennen huoltoa ravistin ja apuvälineet tulisi puhdistaa ja dekontaminoida läpikotaisin jotta täysi ja turvallinen tarkastus voidaan suorittaa.

Thermo Fisher Scientific tarjoaa tarkistus- ja huoltosopimuksia tähän työhön. Kaikki tarpeelliset korjaukset tehdään maksutta takuukauden aikana ja maksua vastaan sen jälkeen. Tämä pätee vain, jos ravistimen on huoltanut valtuutettu Thermo Fisher Scientificin huoltoteknikko.

### 4.10. Toimitus ja hävittäminen

VAROITUS Vaara terveydelle tartuntavaarallisista aineista. Kun ravistin ja apuvälineet poistetaan käytöstä poisheittoa varten, ne täytyy puhdistaa ja jos tarpeen desinfioida ja dekontaminoida. Ongelmatapauksissa ota yhteysThermo Fisher Scientificin asiakaspalveluun.

Ravistimen poisheitossa huomioi maan määräykset. Ota yhteys Thermo Fisher Scientificin asiakaspalveluun ravistimen poisheittoa varten. Yhteystiedot: tämän käyttöohjeen takasivu tai <u>www.thermofisher.com</u>.

Ota huomioon kuljetus- ja toimitustiedot kohdassa ("Kuljetus" sivulla 52 ja kohdassa "Toimitus" sivulla 59).

# 5. Vianmääritys

Ei.	Kuvaus	Ratkaisut	Kuvake
Virheet			
11999	Tapahtui sisäinen virhe.	Tapahtui sisäinen virhe. Käynnistä laite uudelleen katkaisemalla virta ja kytkemällä se takaisin päälle. Jos vikailmoitus yhä näkyy, ota yhteys huoltoteknikkoon.	
Varoitukset			
3	Lämpötilan kalibroinnin määräaika. Se on ollut 1 vuosi.	Yhden vuoden voimassaoloaika on päättynyt. Uudista lämpötilan kalibrointi, ks. osio "4. 6. Lämpötilan kalibrointi" sivulla 126.	CAL
4	Virta on uudelleen päällä ja automaattinen uudelleenkäynnistys mahdollistettu.	Sähkökatkos tapahtui edellisen ajon aikana. Kun virta on saatu takaisin, ajo on automaattisesti käynnistetty uudelleen.	Ę
6	Ympäristön lämpötila-anturit vialliset.	Ympäristön lämpötila-anturin mittaama lämpötila ei ole uskottava.	
10	Puhaltimen 1 nopeus yli rajan.	Puhaltimen 1 nopeus ylittää uskottavuuskynnyksen.	•2
11	Puhaltimen 2 nopeus yli rajan.	Puhaltimen 2 nopeus ylittää uskottavuuskynnyksen.	•2
12	Puhaltimen 3 nopeus yli rajan.	Puhaltimen 3 nopeus ylittää uskottavuuskynnyksen.	•2
13	Puhaltimen 4 nopeus yli rajan.	Puhaltimen 4 nopeus ylittää uskottavuuskynnyksen.	•2
20	Kansi on ollut kauan auki.	Sulje ravistimen kansi, jotta kammion lämpötila säilyy.	
Hälytykset			
1	Korkean lämpötilan hälytys.	Kammion lämpötila on ylittänyt hälytyksen ylärajan. Tarkista näytteet, ympäristöolosuhteet ja/ tai muuta asetuksia.	
2	Matalan lämpötilan hälytys.	Kammion lämpötila on ylittänyt hälytyksen alarajan. Tarkista näytteet, ympäristöolosuhteet ja/tai muuta asetuksia.	
3	Automaattinen uudelleenkäynnistys epäonnistuneen virtakatkon jälkeen.	Sähkökatkos tapahtui edellisen ajon aikana. Automaattista uudelleenkäynnistämistä ei voitu suorittaa onnistuneesti.	Ç,
5	Ajon kiihtyvyys liian hidas. Haluttua nopeutta ei voida saavuttaa.	Haluttua nopeutta ei voinut saavuttaa ajoissa. Tarkista asetuksesi ja/tai lasti (apuvälineet ja näytteet) alustalla.	$\odot$

#### Toimitus ja hävittäminen

Ei.	Kuvaus	Ratkaisut	Kuvake
7	Puhaltimen 1 nopeusmittaus ilmoittaa odottamattomasta pysähtymisestä.	Käännä laite pois päältä ja taas takaisin päälle. Jos virheilmoitus ei poistu, ota yhteys huoltoteknikkoon.	•
8	Puhaltimen 2 nopeusmittaus ilmoittaa odottamattomasta pysähtymisestä.	Käännä laite pois päältä ja taas takaisin päälle. Jos virheilmoitus ei poistu, ota yhteys huoltoteknikkoon.	•
9	Puhaltimen 3 nopeusmittaus ilmoittaa odottamattomasta pysähtymisestä.	Käännä laite pois päältä ja taas takaisin päälle. Jos virheilmoitus ei poistu, ota yhteys huoltoteknikkoon.	•
10	Puhaltimen 4 nopeusmittaus ilmoittaa odottamattomasta pysähtymisestä.	Käännä laite pois päältä ja taas takaisin päälle. Jos virheilmoitus ei poistu, ota yhteys huoltoteknikkoon.	•
12	Ajon käynnistysvirhe – ei saada nopeussignaalia.	Ravistimen alusta on jumissa. Tarkista että alustan ympärillä on tarpeeksi tilaa ja/tai vähennä lastia (apuvälineet ja näytteet) alustalla. Sen jälkeen paina ravistimessa uudestaan START. Jos virheilmoitus ei poistu, ota yhteys huoltoteknikkoon.	
24	Nopeuden mittauksen virhe ajon aikana.	Epänormaali nopeudenmuutos havaittu. Tarkista lasti (apuvälineet ja näytteet) ja/tai puristimet ravistimen alustassa. Sen jälkeen paina ravistimessa uudestaan START. Jos virheilmoitus ei poistu, ota yhteys huoltoteknikkoon.	
26	Nopeusmittaus havaitsi odottamattoman pysäyksen ajon aikana.	Tarkista että alustan ympärillä on tarpeeksi tilaa ja/tai vähennä lastia (apuvälineet ja näytteet) alustalla. Varmista, ettei sulake ravistimen takana ole lauennut ("Sulake" sivulla 47). Sen jälkeen paina ravistimessa uudestaan START. Jos virheilmoitus ei poistu, ota yhteys huoltoteknikkoon.	٢
82	Moottorin virtalukemat osoittavat ylikuormitusta.	Moottorin ylikuormitus havaittu. Älä lastaa tai pura alustaa ajon aikana. Alenna nopeutta tai säädä lastia (apuvälineet ja näytteet) alustalla.	
83	Moottorin virtamittaus ei ole rajoissa.	Moottorin ylikuormitus havaittu. Älä lastaa tai pura alustaa ajon aikana. Alenna nopeutta tai säädä lastia (apuvälineet ja näytteet) alustalla.	

Taulukko 59: Luettelo virheistä, hälytysilmoituksista ja hälytyksistä

HUOMIO Jos näytössä näkyy virheviesti, jota ei ole tässä taulukossa, ota yhteys huoltoteknikkoon.

# GPL (yleinen julkinen lisenssi)

Laitteen ohjelmiston jotkin osat käyttävät avoimen lähdekoodin ohjelmistoja, jotka on julkaistu GPL-, LGPL- tai muun avoimen lähdekoodin lisenssin nojalla. Nämä ovat taulukossa yksityiskohtaisesti luetellut kirjastot. Jos kyseinen lisenssi antaa lähdekoodin käytetyille kirjastoille (kolmansien osapuolten kirjastoille), se on saatavana Thermo Fisher Scientificilta. Käytettävän avoimen lähdekoodin ohjelmiston vastaavat käyttöoikeusehdot ovat osa mukana toimitettua lähdekoodipakettia.

Kirjasto	Versio	Käyttöönotto	Lisenssi
Qt	5.8	BSP	LGPLv3
Log4Cplus	1.2.0	Sovellus	ApachePublicLicense v2/ Twoclause BSD license
boost	1.72.0	Sovellus	Boost License 1.0
json11	1.0.0	Sovellus	MIT License

#### Kolmansien osapuolten kirjastot

\_\_\_\_\_

https://www.gnu.org/licenses/lgpl-3.0.de.html https://www.gnu.org/licenses/lgpl-3.0.en.html

## Sisältö

#### Α

Aika 75 Aikamoodi 73 Alue 76 Alustan asennus 89 Alustat 18 Apuvälineiden tarkistus 123 Asetukset 68 Autoklavointi 125 Automaattinen uudelleenkäynnistys 73 Äänenvoimakkuus 69

### D

Dekantterilasitelineen asennus 101 Dekontaminaatio 124 Desinfiointi 124 Direktiivit 43 Dual Stack-alustat 91

### Ε

Erotussuppilon pidikkeen asennus 102 Ethernet 47

### G

GPL (yleinen julkinen lisenssi) 137 Graafinen käyttöliittymä 60

### Η

Hallintalaitteet 73 Huolto 121 Hälytykset 69 Hävittäminen 134

### K

Kaasunjakoputkiston asennus 113 Kaaviot 86 Kalibraatio 72 Kallistuskulmaltaan säädettävän koeputkitelineen pidikkeen asennus 97 Kieli 74 Kirkkaus 73 Koepullopidikkeen ja astian asennus 92 Koeputkitelineen asennus 96 Korkein nopeus 118 Kuljetus 52 Kunnossapito 77, 121, 134 Käynnistys 54 Käyttö 60 Käyttöohjeessa käytetyt symbolit 7 Käyttötarkoitus 5

### L

Laitteen nimi 76 Lastaus 115 Lataa 116 Liimakiinnitteisen maton asennus ja käyttö 106 Liitännät 46 Lämpötilan asetusarvo 64, 65, 73 Lämpötilan kalibrointi 126 Lämpötilan määritys 64

### Μ

Metalliosat 123 Mikrolevyn / Deepwell-levyn pidikkeen asennus 98 Mukauta valikkoa 76 Muoviosat 123

### Ν

Nelikulmaisen pidikkeen asennus 94 Nopeuden asetuspiste 73 Normaali käyttö 115, 117 Nukkumismoodi 75 Näyttö 73 0

Ohjelmat 78

#### Ρ

Puhdistus 123 Päiväyksen muoto 75 Päiväys 74 Pääsyn valvonta 71

### S

Signaalisanat ja symbolit 6 Sijoituspaikka 50 Standardit 43 Sulake 47 Säädettävän astianpidikkeen asennus 99

### Т

Tapahtumaloki 84 Tarviketason asennus 105 Tarvikkeet 17, 88 Tekniset tiedot 11 Tiedostot ja info 77 Toimitus 59, 134 Torkun aikakatkaisu 70 Tuotekatsaus 44 Tuotteet 49 Turvallinen nopeus 117

### U

**USB** 48

### V

Varoitukset 69 Verkkovirtaliitäntä 47 Vianmääritys 135

#### Υ

Yleisalustat 90

# thermo scientific





Valmistettu

Thermo Electron LED GmbH:lle Zweigniederlassung Osterode Am Kalkberg, 37520 Osterode am Harz Germany Alkuperämaa: USA

CE

Thermo Scientific Solaris 2000 Thermo Scientific Solaris 2000 I Thermo Scientific Solaris 2000 R Thermo Scientific Solaris 4000 Thermo Scientific Solaris 4000 I Thermo Scientific Solaris 4000 R

70900190 on alkuperäinen käyttäjän opas. Tämä käyttöopas on alkuperäisen käyttöoppaan käännös.

#### thermofisher.com/shaker

© 2019–2024 Thermo Fisher Scientific Inc. Kaikki oikeudet pidätetään.

Kaikki tuotenimet ovat Thermo Fisher Scientific Inc.:in ja sen tytäryhtiöiden omistuksessa jos ei toisin mainita. Kaikki tuotteet eivät ole saatavilla kaikissa maissa. Konsultoikaa yksityiskohdista paikallista myyntiedustajaanne.

Näytetyt kuvat käsikirjassa ovat esimerkkejä ja ne voivat erota asetettujen parametrien ja kielen osalta.

Australia +61 39757 4300

**Itävalta** +43 1 801 40 0

Belgia +32 53 73 42 41

**Kiina** +800 810 5118 **tai** +400 650 5118

**Ranska** +33 2 2803 2180

Saksa paikallinen maksuton 0800 1 536 376

**Saksa kansainvälinen** +49 6184 90 6000 Intia +91 22 6716 2200

**Italia** +39 02 95059 552

**Japani** +81 3 5826 1616

Alankomaat +31 76 579 55 55

**Uusi Seelanti** +64 9 980 6700

Pohjoimaat/Baltia/IVY-maat +358 10 329 2200

Veäjä +7 812 703 42 15 Espanja/Portugali +34 93 223 09 18

**Sveitsi** +41 44 454 12 12

Yhdistynyt Kuningaskunta/ Irlanti +44 870 609 9203

**USA/Kanada** +1 866 984 3766

Muut Aasian maat +852 2885 4613

**Listaamattomat maat** +49 6184 90 6000

