

Thermo Scientific rysteapparat i Solaris-serien

Brugsanvisning

70900198-d • 2023-03



Indhold

Forord	
Om denne	e brugsanvisning 5
Reglemer	nteret anvendelse 5
Ū	Åbne rysteapparater 5
	Inkuberede rysteapparater (kun opvarmede) og kølederysteapparater (kølede og opvarmede)5Påkrævet viden5
Signalord	og symboler 6
	Anvendte symboler på rysteapparatet og tilbehørsdelene 6
	Symboler, der anvendes i brugsanvisning 7
Sikkerhed	Isanvisninger7
1. Teknisk	e specifikationer
1.1. Tek	niske data
	Rysteapparat Thermo Scientific Solaris 200011
	Rysteapparat Thermo Scientific Solaris 400012
	Rysteapparat Thermo Scientific Solaris 2000 I13
	Rysteapparat Thermo Scientific Solaris 4000 I14
	Rysteapparat Thermo Scientific Solaris 2000 R15
	Rysteapparat Thermo Scientific Solaris 4000 R16
1. 2. Tilb	ehør
1. 2. 1.	Platforme
1. 2. 2.	Kolbeholder
1. 2. 3.	Kolbeholder til mikrotiter-/Deepwell-plader 23
1. 2. 4.	Rørstativ
1. 2. 5.	Rørstativholder med justerbar vinkel 28
1. 2. 6.	Justerbare beholderholdere 30
1. 2. 7.	Holder til firkantede beholdere
1. 2. 8.	Nalgene bægerglasstativ 34
1. 2. 9.	Skilletragtholder 38
1. 2. 10.	Universalbakke 39
1. 2. 11.	Klæbemåtter 40
1. 2. 12.	Generelt tilbehør 42

1. 3.	Star	ndarder og direktiver	43
1. 4.	Proc	duktoversigt	44
1.	4. 1.	Solaris 2000	44
1.	4. 2.	Solaris 4000	44
1.	4. 3.	Solaris 2000 I / 2000 R	45
1.	4.4.	Solaris 4000 I / 4000 R	45
1.	4. 5.	Tilslutninger	46
2. Trar	nspo	rt og opstilling	49
2. 1.	Udp	akning	49
	•	Leveringsomfang	49
2. 2.	Ops	tillingssted	50
2. 3.	Trar	nsport	52
		Håndtering af rysteapparatet	52
2. 4.	Nive	ellering	53
2. 5.	Lysr	nettilslutning	53
2. 6.	Gru	ndindstilling	54
2. 7.	Opb	evaring	60
2. 8.	Fors	sendelse	60
3. Beti	ienin	a	61
3 1	Tæn	od/sluk	61
2.0			61
J. Z.	01a 2 1	Indetilling of nysteenparatets vigtigste driftsparametre	62
о. З	2. 1. 2. 2	Driftstilstand	67
3.	2.3.	Indstillinger	69
3.	2.4.	Programmer	79
3. 3.	Tilb	- ehør	89
3.	3. 1.	Montering af platformen	90
3.	3. 2.	Isætning af kolbeholder og beholder	93
3.	3. 3.	Installation af firkantet kolbeholder	95
3.	3. 4.	Installation af rørstativ	97
3.	3. 5.	Installation af rørstativholder med justerbar vinkel	98
3.	3. 6.	Installation af kolbeholder til mikrotiter-/Deepwell-plader	99
3.	3. 7.	Installation af justerbar beholderholder	100

3. 3. 8. Installation af bægerglasstativer	102
3. 3. 9. Installation af skilletragtholder	103
3. 3. 10. Installation af universalbakke	106
3. 3. 11. Lægning og anvendelse af klæbemåtten	107
3. 3. 12. Installation af gasfordeler	114
3. 4. Fyldning og tilsigtet anvendelse	116
Påfyldning	117
Reglementeret anvendelse	118
4. Vedligeholdelse og pleje	122
4. 1. Principper	123
Kontrol af tilbehøret	124
4. 2. Rengøring	124
Kontrolpanel	124
4. 3. Desinficering	125
4. 4. Dekontaminering	125
4. 5. Autoklavering	126
4. 6. Temperaturkalibrering	127
4. 6. 1. Pålidelige metoder til temperaturkalibrering	127
4. 6. 2. Gennemførelse af temperaturkalibrering	127
4. 6. 3. Temperaturkalibreringsforespørgsel	131
4.7. Firmwareinstallation	131
4.8. Udskiftning af platformen	135
4. 9. Service	135
1 10 Earaandalaa ay bartakaffalaa	135
5. Årsager til fejl og fejlafhjælpning	136
5. Årsager til fejl og fejlafhjælpning	136 138
5. Årsager til fejl og fejlafhjælpning GPL (General Public License, offentlig licens)	136 138 139

Forord

Om denne brugsanvisning

Denne brugsanvisning beskriver følgende rysteapparater og de tilhørende tilbehørsdele ("1. 2. Tilbehør" på side 17):

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

Beskrivelserne og anvisningerne i denne brugsanvisning relaterer ikke til det fuldstændige produktnavn, men til de enkelte rysteapparater efter klasser:

- Åbne rysteapparater
- Inkuberede rysteapparater (kun opvarmede)
- Kølede rysteapparater (kølede og opvarmede)

Reglementeret anvendelse

Åbne rysteapparater

Dette apparat anvendes inden for forskning og fremstilling til at ryste opløsninger i en cirkulær 2D-bane og samtidig styre anvendelsens omdrejningstal og varighed.

Inkuberede rysteapparater (kun opvarmede) og kølede

rysteapparater (kølede og opvarmede)

Dette apparat anvendes inden for forskning og fremstilling til at ryste opløsninger i en cirkulær 2D-bane og samtidig styre anvendelsens omdrejningstal, varighed og temperatur.

Påkrævet viden

Dette rysteapparat må kun betjenes af uddannet personale.

Signalord og symboler

Signalord	Risikoniveau
ADVARSEL	Henviser til farlige situationer, der kan medføre dødbringende eller alvorlige kvæstelser, hvis de ikke undgås.
FORSIGTIG	Henviser til farlige situationer, der kan medføre lettere eller moderate kvæstelser, hvis de ikke undgås.
BEMÆRK	Henviser til vigtige informationer, der ikke er forbundet med farer.

Anvendte symboler på rysteapparatet og tilbehørsdelene

Løs og følg anvisningerne i brugsanvisningen, så du ikke udsætter dig selv eller dine omgivelser for fare.

8	Du finder yderligere oplysninger om dette i brugsanvisningen
	Træk netstikket ud
	Generel risiko
	Risiko for forbrænding på grund af varme overflader
	Fare for finger- eller håndskader på grund af mekaniske dele, der lukker utilsigtet

Symboler, der anvendes i brugsanvisning

Løs og følg anvisningerne i brugsanvisningen, så du ikke udsætter dig selv eller dine omgivelser for fare.

Generel risiko	4	Risiko for elektrisk stød
Biologisk betinget fare	\bigwedge	Fare for snitsår
Risiko på grund af brændbare materialer	í	Henviser til vigtige informationer, der ikke er forbundet med farer.
Fare for finger- eller håndskader på grund af mekaniske dele, der lukker utilsigtet.		

Sikkerhedsanvisninger



ADVARSEL

Hvis disse sikkerhedsanvisninger tilsidesættes, kan der opstå farlige situationer, der kan medføre dødbringende eller alvorlige kvæstelser, hvis sådanne situationer ikke undgås.

- Læs og følg sikkerhedsanvisningerne.
- Rysteapparatet må kun anvendes til dets tilsigtede anvendelse. Hvis den ikke anvendes til det tilsigtede formål, kan dette medføre skader, kontaminering og livsfarlige kvæstelser.
- Rysteapparatet må kun betjenes af uddannet personale.
- Det er driftsherrens pligt at sikre, at der anvendes egnede personlige værnemidler. Læs og følg Verdenssundhedsorganisationen WHO's Laboratory Biosafety Manual og nationale bestemmelser.



Skader på grund af forkert strømforsyning.

Sørg for, at rysteapparatet kun tilsluttes til korrekt jordforbundne stikkontakter.

ADVARSEL



Fare ved håndtering af farlige stoffer.

ADVARSEL

Hvis du arbejder med korrosive prøver (saltopløsninger, syrer, baser), skal du rengøre tilbehørsdelene og rysteapparatet omhyggeligt.

- Rysteapparatet er hverken inertiseret eller eksplosionsbeskyttet. Brug aldrig rysteapparatet i eksplosionsfarlige omgivelser.
- Ryst ikke giftige eller radioaktive materialer samt patogene mikroorganismer uden passende sikkerhedsforanstaltninger.
- Læs og følg Verdenssundhedsorganisationen WHO's Laboratory Biosafety Manual og nationale bestemmelserne, hvis du ryster nogen form for farlige materialer. Hvis der rystes mikrobiologiske prøver af risikogruppe II (iht. Verdenssundhedsorganisationen WHO's Laboratory Biosafety Manual), skal der anvendes aerosoltætte biotætninger. Du finder "Laboratory Biosafety Manual" på Verdenssundhedsorganisationens internetside (www.who.int). For materialer i en højere risikogruppe skal der træffes yderligere sikkerhedsforanstaltninger.
- Hvis giftige eller patogene substanser har kontamineret rysteapparatet eller dele af det, skal du gennemføre passende desinfektions- og dekontamineringsforanstaltninger ("Dekontaminering" på side 125; "Desinficering" på side 125).
- Afbryd strømforsyningen til rysteapparatet og forlad omgående området, hvis der opstår en farlig situation.



ADVARSEL

Sundhedsskader på grund af smitsomme stoffer.

Hvis der utilsigtet kommer væske eller andre materialer under platformen, skal rysteapparatet omgående slukkes, lysnetstikket skal trækkes ud, og platformen skal fjernes ("Platforme" på side 18).

Tør spildt væske op iht. de gældende standardprocedurer for laboratorierne. Anvend korrekte personlige værnemidler.



Sundhedsskader på grund af at der rystes eksplosive eller brændbare materialer eller stoffer.

ADVARSEL

Ryst ikke eksplosive eller brændbare materialer eller stoffer.



Forbrænding på varme overflader.

FORSIGTIG

I et temperaturstyret rysteapparat kan overflader som platformen og tilbehørsdele under låget blive meget varme. Rør ikke ved varme overflader. Vent, indtil de varme overflader er afkølet.



FORSIGTIG

Snitsår på grund af glasskår.

En løsnet platform og tilbehørsdele, som f.eks. kolbeholderen, kan medføre, at glasset går i stykker, hvis beholderne falder ud af rysteapparatet. Sørg for, at platformen og tilbehørsdelene installeres korrekt ved at anvende de korrekte værktøjer og skruer. Kontroller, at kolbeholderne er designet til at holde beholdernes og prøvernes vægt ved det valgte omdrejningstal. Kolbeholdere med individuel påfyldning er muligvis kun egnet til omdrejningstal under rysteapparatets stabilitetsgrænseværdier. Du finder yderligere oplysninger i afsnittet "3. 4. Fyldning og tilsigtet anvendelse" på side 116. Vær opmærksom på usædvanlige lyde. De kan stamme fra en løsnet platform og tilbehørsdele.



Beskadigelse af rysteapparatet eller funktionsfejl på grund af et beskadiget betjeningspanel.

- Brug ikke rysteapparatet, hvis betjeningspanelet er beskadiget.
- Sluk rysteapparatet. Træk netstikket ud. Lad en autoriseret servicetekniker udskifte betjeningspanelet.



FORSIGTIG

Biologisk risiko på grund af ødelagte eller utætte beholdere.

Ukorrekt installerede tilbehørsdele kan medføre, at prøverne bliver spildt.

- Kontroller, at tilbehørsdelene er installeret korrekt ved at anvende de rigtige værktøjer og skruer.
- Kontroller, at tilbehørsdelene er placeret korrekt på platformen.
- Anvend altid kun beholdere med tilbehør, der er tilpasset den pågældende størrelse.
- Beholderne skal være intakte og installeret korrekt.

FORSIGTIG



Sikkerheden kan blive forringet på grund af forkert fyldning og beskadigede tilbehørsdele.

FORSIGTIG

- Sørg altid for en så ensartet lastfordeling som mulig (tilbehør og prøver), især ved brug af en totrins platform.
- Brug aldrig tilbehørsdele, der viser tegn på korrosion, eller som har revner. Kontakt kundeservice for yderligere oplysninger.
- Anvend kun rysteapparatet, når det er fyldt korrekt.
- Overfyld aldrig rysteapparatet.
- Kontroller, at tilbehøret er installeret korrekt, før du tager rysteapparatet i brug. Følg anvisningerne i afsnittet "Tilbehør" på side 89.



FORSIGTIG

Risiko for personskade ved manglende overholdelse af funktionsprincipperne.

- Anvend aldrig rysteapparatet uden en korrekt monteret platform.
- Anvend aldrig rysteapparatet, hvis dele af beklædningen er beskadiget eller fjernet.
- Anvend aldrig et rysteapparat med en defekt lågfjeder.
 En defekt lågfjeder kan ikke holde låget sikkert i helt åben position.
- Bevæg ikke rysteapparatet mens det er i drift.
- Læn dig ikke op ad rysteapparatet.
- Fyld eller tøm altid først rysteapparatet når det står helt stille, og dette vises på betjeningspanelet.
- Læg ikke noget på rysteapparatet mens det er i drift.
- Rør ikke ved platformen eller tilbehørsdelene på rysteapparatet mens det er i drift.
- Rysteapparatets hus må ikke åbnes af brugeren.



Beskyttelsen kan forringes ved brug af ikke-kompatible tilbehørsdele.

BEMÆRK

Anvend kun tilbehørsdele til dette rysteapparat, der er godkendt af Thermo Fisher Scientific. Du finder de opdaterede lister på <u>www.thermofisher.com</u>.



Sådan slukkes rysteapparatet:

Tryk på knappen STOP. Sluk rysteapparatet med hovedafbryderen. Træk netstikket ud. Afbryd strømtilførslen i en nødsituation.

BEMÆRK

1. Tekniske specifikationer

1.1. Tekniske data

Rysteapparat Thermo Scientific Solaris 2000					
Omdrejningstalområde	15 - 525 o/min				
Driftstid	99 timer 59 min (i trin på	1 min) eller kontinuerlig dr	ift		
Støjniveau ved maks. omdrejningstal	50 dB (A) (1 m foran app	paratet i en højde på 1,6 m)		
Maks. fyldning (inkl. platform, tilbehør og prøver)	25 kg (55 lbs)	25 kg (55 lbs)			
Elektrisk tilslutning	100–240 V ±10 %, 50 /	60 Hz			
Strømforbrug	60 W				
Omgivelsesbetingelser					
Ved opbevaring og forsendelse	Temperatur: -10 °C til 58 Fugtighed: 15 % til 85 %	5 °C			
l drift	Andendelse i indendørs rum Højde op til 3 000 m over normal nul Temperatur: 5 °C til 40 °C Maks. relativ fugtighed 80 % til 31 °C, lineært aftagende til 50 % relativ fugtighed ved 40 °C.				
Forureningsgrad	2				
Overspændingskategori	I				
IP	20				
Grænseflader	USB	Ethernet			
	2x USB-A 2.0	RJ45			
Dimensioner	Længde	Bredde	Højde		

	47 cm (18,5 in)	37 cm (14,5 in)	15 cm (5,5 in)
Vast	20.0 kg (46.0 lbs)		
Vægt	20,8 Kg (40,0 108)		

Tabel 1: Tekniske data Solaris 2000

Rysteapparat Thermo Scientific Solaris 4000

Omdrejningstalområde	15 - 525 o/min			
Driftstid	99 timer 59 min (i trin på 1 min) eller kontinuerlig drift			
Støjniveau ved maks. omdrejningstal	54 dB (A) (1 m foran apparatet i en højde på 1,6 m)			
Maks. fyldning (inkl. platform, tilbehør og prøver)	43 kg (95 lbs)			
Elektrisk tilslutning	100–240 V ±10 %, 50 /	60 Hz		
Strømforbrug	80 V			
Omgivelsesbetingelser				
Ved opbevaring og forsendelse	Temperatur: -10 °C til 55 Fugtighed: 15 % til 85 %	5 °C		
l drift	Andendelse i indendørs rum Højde op til 3 000 m over normal nul Temperatur: 5 °C til 40 °C Maks. relativ fugtighed 80 % til 31 °C, lineært aftagende til 50 % relativ fugtighed ved 40 °C.			
Forureningsgrad	2			
Overspændingskategori	11			
IP	20			
Grænseflader	USB	Ethernet		
	2x USB-A 2.0	RJ45		
Dimensioner	Længde	Bredde	Højde	
	65 cm (25,6 in)	58 cm (22,8 in)	18 cm (7 in)	
Vægt	75,1 kg (165,5 lbs)			

Tabel 2: Tekniske data Solaris 4000

Rysteapparat Thermo Scientific Solaris 2000 I

Omdrejningstalområde	15 - 525 o/min		
Temperaturindstillingsområde	30-60 °C		
Temperaturstabilitet i kolben ved 37 °C	Omgivelsestemperatur 23 °C i 1 time ved en stabil temperatur i prøverummet \pm 0,1 °C		
Temperaturensartethed i kolben ved 37 $^{\circ}\mathrm{C}$	Omgivelsestemperatur 23 ± 0,5 °C	°C i 1 time ved en stabil tem	peratur i prøverummet
Driftstid	99 timer 59 min (i trin på	1 min) eller kontinuerlig dri	ift
Støjniveau ved maks. omdrejningstal	52 dB (A) (1 m foran app	aratet i en højde på 1,6 m)
Maks. fyldning (inkl. platform, tilbehør og prøver)	16 kg (35 lbs), inkl. 11 x 14 platform på	2,3 kg (5,1 lbs)	
Elektrisk tilslutning	100–120, 200–240 V ±	10 %, 50 / 60 Hz	
Strømforbrug	900 W		
Omgivelsesbetingelser			
Ved opbevaring og forsendelse	Temperatur: -10 °C til 55 Fugtighed: 15 % til 85 %	°C	
l drift	Andendelse i indendørs rum Højde op til 3 000 m over normal nul Temperatur: 5 °C til 40 °C Maks. relativ fugtighed 80 % til 31 °C, lineært aftagende til 50 % relativ fugtighed ved 40 °C.		
Forureningsgrad	2		
Overspændingskategori			
P	20		
Grænseflader	USB	Ethernet	
	2x USB-A 2.0	RJ45	
Dimensioner	Længde	Bredde	Højde
	70 cm (27,6 in)	36 cm (14,2 in)	46 cm (18,1 in)
Med åbent låg			79 cm (30,8 in)
Vægt	44,9 kg (99 lbs)		

Tabel 3: Tekniske data Solaris 2000 I

Rysteapparat Thermo Scientific Solaris 4000 I

Omdrejningstalområde	15 - 525 o/min		
Temperaturindstillingsområde	30–60 °C		
Temperaturstabilitet i kolben ved 37 °C	Omgivelsestemperatur 23 °C i 1 time ved en stabil temperatur i prøverummet \pm 0,1 °C		
Temperaturensartethed i kolben ved 37 °C	Omgivelsestemperatur 23 ' ± 0,5 °C	°C i 1 time ved en stabil temp	ceratur i prøverummet
Driftstid	99 timer 59 min (i trin på	1 min) eller kontinuerlig dri	ft
Støjniveau ved maks. omdrejningstal	52 dB (A) (1 m foran app	aratet i en højde på 1,6 m)
Maks. fyldning (inkl. platform, tilbehør og prøver)	32 kg (70,5 lbs), inkl. 18 x 18 platform på	4,3 kg (9,5 lbs)	
Elektrisk tilslutning	100–120, 200–240 V ±1	0 %, 50 / 60 Hz	
Strømforbrug	900 W		
Omgivelsesbetingelser			
Ved opbevaring og forsendelse	Temperatur: -10 °C til 55 Fugtighed: 15 % til 85 %	°C	
l drift	Andendelse i indendørs rum Højde op til 3 000 m over normal nul Temperatur: 5 °C til 40 °C Maks. relativ fugtighed 80 % til 31 °C, lineært aftagende til 50 % relativ fugtighed ved 40 °C.		
Forureningsgrad	2		
Overspændingskategori	1		
P	20		
Grænseflader	USB	Ethernet	
	2x USB-A 2.0	RJ45	
Dimensioner	Længde	Bredde	Højde
	77 cm (30,3 in)	57 cm (22,4 in)	55 cm (21,7 in)
Med åbent låg			94 cm (36,7 in)
Vægt	68,1 kg (150,1 lbs)		

Tabel 4: Tekniske data Solaris 4000 I

Rysteapparat Thermo Scientific Solaris 2000 R

Omdrejningstalområde	15 - 525 o/min				
Temperaturindstillingsområde	5–60 °C	5–60 °C			
Temperaturstabilitet i kolben ved 37 °C	Omgivelsestemperatur 23 °C i 1 time ved en stabil temperatur i prøverummet \pm 0,1 °C				
Temperaturensartethed i kolben ved 37 $^{\circ}\mathrm{C}$	Omgivelsestemperatur 23 ± 0,5 °C	°C i 1 time ved en stabil tem	peratur i prøverummet		
Driftstid	99 timer 59 min (i trin på	1 min) eller kontinuerlig dri	ift		
Støjniveau ved maks. omdrejningstal	56 dB (A) (1 m foran app	aratet i en højde på 1,6 m)		
Maks. fyldning (inkl. platform, tilbehør og prøver)	16 kg (35 lbs), inkl. 11 x 14 platform på	2,3 kg (5,1 lbs)			
Elektrisk tilslutning	100–240 V ±10 %, 50 /	60 Hz			
Strømforbrug	350 W				
Omgivelsesbetingelser					
Ved opbevaring og forsendelse	Temperatur: -10 °C til 55 °C Fugtighed: 15 % til 85 %				
l drift	Andendelse i indendørs rum Højde op til 3 000 m over normal nul Temperatur: 5 °C til 40 °C Maks. relativ fugtighed 80 % til 31 °C, lineært aftagende til 50 % relativ fugtighed ved 40 °C.				
Forureningsgrad	2				
Overspændingskategori	1				
P	20				
Grænseflader	USB	Ethernet			
	2x USB-A 2.0	RJ45			
Dimensioner	Længde	Bredde	Højde		
	70 cm (27,6 in)	36 cm (14,2 in)	46 cm (18,1 in)		
Med åbent låg			79 cm (30,8 in)		
Vægt	47,2 kg (104,1 lbs)				

Tabel 5: Tekniske data Solaris 2000 R

Rysteapparat Thermo Scientific Solaris 4000 R

Omdrejningstalområde	15 - 525 o/min			
Temperaturindstillingsområde	4–60 °C			
Temperaturstabilitet i kolben ved 37 °C	Omgivelsestemperatur 23 °C i 1 time ved en stabil temperatur i prøverummet \pm 0,1 °C			
Temperaturensartethed i kolben ved 37 °C	Omgivelsestemperatur 23 % ± 0,5 °C	C i 1 time ved en stabil temp	eratur i prøverummet	
Driftstid	99 timer 59 min (i trin på 1	min) eller kontinuerlig driff	t	
Støjniveau ved maks. omdrejningstal	56 dB (A) (1 m foran appa	rratet i en højde på 1,6 m)		
Maks. fyldning (inkl. platform, tilbehør og prøver)	32 kg (70,5 lbs), inkl. 18 x 18 platform på 4	I,3 kg (9,5 lbs)		
Elektrisk tilslutning	100–240 V ±10 %, 50 / 6	60 Hz		
Strømforbrug	750 W			
Omgivelsesbetingelser				
Ved opbevaring og forsendelse	Temperatur: -10 °C til 55 °C Fugtighed: 15 % til 85 %			
l dríft	Andendelse i indendørs rum Højde op til 3 000 m over normal nul Temperatur: 5 °C til 40 °C Maks. relativ fugtighed 80 % til 31 °C, lineært aftagende til 50 % relativ fugtighed ved 40 °C.			
Forureningsgrad	2			
Overspændingskategori	1			
IP	20			
Grænseflader	USB	Ethernet		
	2x USB-A 2.0	RJ45		
Dimensioner	Længde	Bredde	Højde	
	77 cm (30,3 in)	57 cm (22,4 in)	55 cm (21,7 in)	
Med åbent låg			94 cm (36,7 in)	
Vægt	74,9 kg (165,1 lbs)			

Tabel 6: Tekniske data Solaris 4000 R

1.2. Tilbehør



Beskyttelsen kan forringes ved brug af ikke-kompatible tilbehørsdele.

FORSIGTIG Anvend kun tilbehørsdele til dette rysteapparat, der er godkendt af Thermo Fisher Scientific.

Du finder de opdaterede lister på www.thermofisher.com.

1.2.1. Platforme

Platform		
Solaris 2000	Vægt	
Universalplatform Thermo Scientific Solaris 12x14	2,7 kg (5,9 lbs)	SK1214
Totrins universalplatform Thermo Scientific Solaris 12x14	6,4 kg (14,1 lbs)	SK1214D
Opgraderingssæt til totrins universalplatform Thermo Scientific Solaris 12x14		SK1214DK
Universalplatform Thermo Scientific Solaris 18x18	4,3 kg (9,4 lbs)	SK1818
Totrins universalplatform Thermo Scientific Solaris 18x18	9,7 kg (21,3 lbs)	SK1818D
Opgraderingssæt til totrins universalplatform Thermo Scientific Solaris 18x18		SK1818DK
Universalplatform Thermo Scientific Solaris 18x24	6,0 kg (13,2 lbs)	SK1824
Solaris 4000		
Universalplatform Thermo Scientific Solaris 18x30	7,5 kg (16,5 lbs)	SK1830
Totrins universalplatform Thermo Scientific Solaris 18x30	16,3 kg (35,9 lbs)	SK1830D
Opgraderingssæt til totrins universalplatform Thermo Scientific Solaris 18x30		SK1830DK
Universalplatform Thermo Scientific Solaris 36x24	12,0 kg (26,4 lbs)	SK3624
Solaris 2000 I / 2000 R		
Universalplatform Thermo Scientific Solaris 11x14	2,3 kg (5,1 lbs)	SK1114
Solaris 4000 I / 4000 R		
Universalplatform Thermo Scientific Solaris 18x18	4,3 kg (9,5 lbs)	SK1818
Eftermonteringssæt og tilbehør		
Kolbeholder-reservedelssæt (skruer)		SK1001
Platform-reservedelssæt, lille (platformsskruer, værktøj)		SK0100
Platform-reservedelssæt, stort (platformsskruer, værktøj)		SK0101
Skruetrækker til tilbehørsdele		75004131

Tabel 7: Tilgængelige platforme

1. 2. 2. Kolbeholder



Til Solaris 2000 platforme

	Artikelnr.	Universal			Artikelnr. Universal Totrin			rins
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18		
10 ml-Erlenmeyerkolbe	30150BI	72	113	157	140	226		
25 ml-Erlenmeyerkolbe	30151	42	64	80	80	124		
50 ml-Erlenmeyerkolbe	30152BI	42	64	80	80	124		
125 ml-Erlenmeyerkolbe	30153	15	32	40	30	52		
250 ml-Erlenmeyerkolbe	30154BI	9	16	24	16	32		
300 ml-Erlenmeyerkolbe	30155	9	16	20	16	32		
500 ml-Erlenmeyerkolbe	30156BI	9	16	20	16	32		
1 I-Erlenmeyerkolbe	30157BI	4	9	10	8	16		
2 I-Erlenmeyerkolbe	30158	3	5	6	-	-		
4 I-Erlenmeyerkolbe	30159	1	4	4	-	-		
5 I-Erlenmeyerkolbe	30159B	1	2	4	-	-		
6 I-Erlenmeyerkolbe	30160	1	2	2	-	-		
2800 ml-Fernbachkolbe	30162	1	4	4	-	-		
2500 ml-kulturkolbe, lav form	30161	1	1	2	-	-		

Tabel 8: Tilgængelige kolbeholdere til Solaris 2000 platforme

Til Solaris 4000 platforme

	Artikelnr.	Universal		Totrins
		18 x 30	36 x 24	18 x 30
10 ml-Erlenmeyerkolbe	30150BI	203	187	402
25 ml-Erlenmeyerkolbe	30151	112	187	220
50 ml-Erlenmeyerkolbe	30152BI	112	187	220
125 ml-Erlenmeyerkolbe	30153	46	83	92
250 ml-Erlenmeyerkolbe	30154BI	28	40	56
300 ml-Erlenmeyerkolbe	30155	28	40	56
500 ml-Erlenmeyerkolbe	30156BI	28	40	56
1 I-Erlenmeyerkolbe	30157BI	14	20	28
2 I-Erlenmeyerkolbe	30158	6	11	12
4 I-Erlenmeyerkolbe	30159	6	8	-
5 I-Erlenmeyerkolbe	30159B	5	8	-
6 I-Erlenmeyerkolbe	30160	3	6	-
2800 ml-Fernbachkolbe	30162	6	8	12
2500 ml-kulturkolbe, lav form	30161	3	6	6

Tabel 9: Tilgængelige kolbeholdere til Solaris 4000 platforme

Til Solaris 2000 I / 2000 R platforme

	Artikelnr.	Universal
		11 x 14
10 ml-Erlenmeyerkolbe	30150BI	59
25 ml-Erlenmeyerkolbe	30151	35
50 ml-Erlenmeyerkolbe	30152BI	35
125 ml-Erlenmeyerkolbe	30153	15
250 ml-Erlenmeyerkolbe	30154BI	7
300 ml-Erlenmeyerkolbe	30155	7
500 ml-Erlenmeyerkolbe	30156BI	7
1 I-Erlenmeyerkolbe	30157BI	4
2 I-Erlenmeyerkolbe	30158	2
4 I-Erlenmeyerkolbe	30159	-
5 I-Erlenmeyerkolbe	30159B	-
6 I-Erlenmeyerkolbe	30160	-
2800 ml-Fernbachkolbe	30162	1
2500 ml-kulturkolbe, lav form	30161	1

Tabel 10: Tilgængelige kolbeholdere til Solaris 2000 I / 2000 R platforme

Til Solaris 4000 I / 4000 R platforme

	Artikelnr.	Universal
		18 x 18
10 ml-Erlenmeyerkolbe	30150BI	113
25 ml-Erlenmeyerkolbe	30151	64
50 ml-Erlenmeyerkolbe	30152BI	64
125 ml-Erlenmeyerkolbe	30153	32
250 ml-Erlenmeyerkolbe	30154BI	16
300 ml-Erlenmeyerkolbe	30155	16
500 ml-Erlenmeyerkolbe	30156BI	16
1 I-Erlenmeyerkolbe	30157BI	9
2 I-Erlenmeyerkolbe	30158	6
4 I-Erlenmeyerkolbe	30159	-
5 I-Erlenmeyerkolbe	30159B	-
6 I-Erlenmeyerkolbe	30160	-
2800 ml-Fernbachkolbe	30162	4
2500 ml-kulturkolbe, lav form	30161	1

Tabel 11: Tilgængelige kolbeholdere til Solaris 4000 I / 4000 R platforme

1. 2. 3. Kolbeholder til mikrotiter-/Deepwell-plader



Til Solaris 2000 platforme

	Artikelnr.	Universal		Totrins		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Mikrotiter- / Deepwell-plade	30175	7	12	16	14	24

Tabel 12: Kolbeholderne til mikrotiter-/Deepwell-pladers kapacitet til Solaris 2000 platforme

Til Solaris 2000 I / 2000 R platforme

	Artikelnr.	Universal
		11 x 14
Mikrotiter- / Deepwell-plade	30175	6

Tabel 13: Mikrotiter-/Deepwell-pladernes kapacitet til Solaris 2000 I / 2000 R platforme

Til Solaris 4000 platforme

	Artikelnr.	Unive	Universal	
		18 x 30	36 x 24	18 x 30
Mikrotiter- / Deepwell-plade	30175	22	36	42

Tabel 14: Mikrotiter-/Deepwell-pladernes kapacitet til Solaris 4000 platforme

Til Solaris 4000 I / 4000 R platforme

	Artikelnr.	Universal
		18 x 18
Mikrotiter- / Deepwell-plade	30175	10

Tabel 15: Mikrotiter-/Deepwell-pladernes kapacitet til Solaris 4000 I / 4000 R platforme

1. 2. 4. Rørstativ

Til Solaris 2000 platforme



	Artikelnr.	Universal		Totrins		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Halv størrelse						
10–13 mm, rød, 6 x 6-anordning	30181	8	12	15	14	24
14–16 mm, orange, 6 x 6-anordning	30183	5	9	11	9	17
17–20 mm, hvid, 4 x 5-anordning	30185	7	11	14	13	21
21–25 mm, blå, 4 x 4-anordning	30187	6	9	11	10	17
26–30 mm, grøn, 3 x 3-anordning	30189	6	9	12	11	18
til Mikrotiterrør, 1,5 ml, blå, 4 x 6-anordning	30191	6	10	13	11	19
Størrelse						
10–13 mm, 6 x 12-anordning	30180BI	3	7	10	6	13
14–16 mm, 6 x 12-anordning	30182	3	4	6	6	8
17–20 mm, 4 x 10-anordning	30184	3	5	7	6	8
21–25 mm, 4 x 10-anordning	30186	2	3	5	4	6
26–30 mm, 3 x 8-anordning	30188	3	4	6	5	8
til Mikrotiterrør, 1,5 ml, 8 x 12-anordning	30190	3	5	7	6	9

Tabel 16: Tilgængelige rørstativer til Solaris 2000 platforme

Til Solaris 4000 platforme

	Artikelnr.	Universal		Totrins	
		18 x 30	36 x 24	18 x 30	
Halv størrelse					
10–13 mm, rød, 6 x 6-anordning	30181	21	32	40	
14–16 mm, orange, 6 x 6-anordning	30183	14	20	26	
17–20 mm, hvid, 4 x 5-anordning	30185	18	20	34	
21–25 mm, blå, 4 x 4-anordning	30187	14	22	25	
26–30 mm, grøn, 3 x 3-anordning	30189	15	24	30	
til Mikrotiterrør, 1,5 ml, blå, 4 x 6-anordning	30191	18	24	34	
Størrelse					
10–13 mm, 6 x 12-anordning	30180BI	14	20	26	
14–16 mm, 6 x 12-anordning	30182	9	12	16	
17–20 mm, 4 x 10-anordning	30184	9	15	18	
21–25 mm, 4 x 10-anordning	30186	7	9	13	
26–30 mm, 3 x 8-anordning	30188	7	10	13	
til Mikrotiterrør, 1,5 ml, 8 x 12-anordning	30190	9	12	17	

Tabel 17: Tilgængelige rørstativer til Solaris 4000 platforme

Til Solaris 2000 I / 2000 R platforme

	Artikelnr.	Universal
Halv størrelse		
10–13 mm, rød, 6 x 6-anordning	30181	6
14–16 mm, orange, 6 x 6-anordning	30183	2
17–20 mm, hvid, 4 x 5-anordning	30185	4
21–25 mm, blå, 4 x 4-anordning	30187	2
26–30 mm, grøn, 3 x 3-anordning	30189	4
til Mikrotiterrør, 1,5 ml, blå, 4 x 6-anordning	30191	3
Størrelse		
10–13 mm, 6 x 12-anordning	30180BI	3
14–16 mm, 6 x 12-anordning	30182	2
17–20 mm, 4 x 10-anordning	30184	2
21–25 mm, 4 x 10-anordning	30186	1
26–30 mm, 3 x 8-anordning	30188	2
til Mikrotiterrør, 1,5 ml, 8 x 12-anordning	30190	2

Tabel 18: Tilgængelige rørstativer til Solaris 2000 I / 2000 R platforme

Til Solaris 4000 I / 4000 R platforme

	Artikelnr.	Universal
		18 x 18
Halv størrelse		
10–13 mm, rød, 6 x 6-anordning	30181	12
14–16 mm, orange, 6 x 6-anordning	30183	8
17–20 mm, hvid, 4 x 5-anordning	30185	10
21–25 mm, blå, 4 x 4-anordning	30187	8
26–30 mm, grøn, 3 x 3-anordning	30189	9
til Mikrotiterrør, 1,5 ml, blå, 4 x 6-anordning	30191	8
Størrelse		
10–13 mm, 6 x 12-anordning	30180BI	7
14–16 mm, 6 x 12-anordning	30182	4
17–20 mm, 4 x 10-anordning	30184	5
21–25 mm, 4 x 10-anordning	30186	3
26–30 mm, 3 x 8-anordning	30188	4
til Mikrotiterrør, 1,5 ml, 8 x 12-anordning	30190	5

Tabel 19: Tilgængelige rørstativer til Solaris 4000 I / 4000 R platforme

1. 2. 5. Rørstativholder med justerbar vinkel

Til Solaris 2000 platforme



	Artikelnr.	Universal			Totrins		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18	
10-13 mm, til 72 prøverør	236090	3	4	6	6	8	
16-20 mm, til 40 prøverør	236091	3	4	6	5	8	
21-25 mm, til 40 prøverør	236092	3	3	4	5	6	
26-30 mm, til 24 prøverør	236093	3	3	4	5	6	

Tabel 20: Tilgængelige rørstativholdere med justerbar vinkel til Solaris 2000 platforme

Til Solaris 4000 platforme

		Artikelnr.	Universal		Totrins	
			18 x 30	36 x 24	18 x 30	
1	0-13 mm, til 72 prøverør	236090	7	14	14	
1	6-20 mm, til 40 prøverør	236091	7	14	14	
2	21-25 mm, til 40 prøverør	236092	5	9	10	
2	26-30 mm, til 24 prøverør	236093	6	9	12	

Tabel 21: Tilgængelige rørstativholdere med justerbar vinkel til Solaris 4000 platforme

Til Solaris 2000 I / 2000 R platforme

	Artikelnr.	Universal
		11 x 14
10-13 mm, til 72 prøverør	236090	2
16-20 mm, til 40 prøverør	236091	2
21-25 mm, til 40 prøverør	236092	1
26-30 mm, til 24 prøverør	236093	1

Tabel 22: Tilgængelige rørstativholdere med justerbar vinkel til Solaris 2000 I / 2000 R platforme

Til Solaris 4000 I / 4000 R platforme

	Artikelnr.	Universal
		18 x 18
10-13 mm, til 72 prøverør	236090	4
16-20 mm, til 40 prøverør	236091	3
21-25 mm, til 40 prøverør	236092	3
26-30 mm, til 24 prøverør	236093	3

Tabel 23: Tilgængelige rørstativholdere med justerbar vinkel til Solaris 4000 I / 4000 R platforme

1. 2. 6. Justerbare beholderholdere





Tabel 24: Tilgængelige justerbare beholderholdere til Solaris 2000 platforme

Til Solaris 4000 platforme

	Artikelnr.	Universal		Totrins	
		18 x 30	36 x 24	18 x 30	
Enkel højdejusterbar beholderholder 11"x14"	75004104	-	-	-	
Dobbelt højdejusterbar beholderholder 12"x14"	75004102	2	2	4	
Dobbelt højdejusterbar beholderholder 18"x18"	75004103	1	2	2	
Justerbar enkelt beholderholder	75004101	8	12	16	

Tabel 25: Justerbare beholderholdere til Solaris 4000 platforme





Til Solaris 2000 I / 2000 R platforme

	Artikelnr.	Universal
		11 x 14
Enkel højdejusterbar beholderholder 11"x14"	75004104	1
Dobbelt højdejusterbar beholderholder 12"x14"	75004102	-
Dobbelt højdejusterbar beholderholder 18"x18"	75004103	-
Justerbar enkelt beholderholder	75004101	2

Tabel 26: Justerbare beholderholdere til Solaris 2000 I / 2000 R platforme

Til Solaris 4000 I / 4000 R platforme

	Artikelnr.	Universal
		18 x 18
Enkel højdejusterbar beholderholder 11"x14"	75004104	-
Dobbelt højdejusterbar beholderholder 12"x14"	75004102	1
Dobbelt højdejusterbar beholderholder 18"x18"	75004103	1
Justerbar enkelt beholderholder	75004101	4

Tabel 27: Justerbare beholderholdere til Solaris 4000 I / 4000 R platforme

1. 2. 7. Holder til firkantede beholdere

Til Solaris 2000 platforme



	Artikelnr.	Universal		Totrins		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Holder til firkantede flasker, 125 ml	75004106	12	16	24	21	32
Holder til firkantede flasker, 250 ml	75004107	9	16	22	18	32
Holder til firkantede flasker, 500 ml	75004108	6	9	12	11	18
Holder til firkantede flasker, 1 000 ml	75004109	4	9	11	8	18

Tabel 28: Tilgængelige holdere til firkantede beholdere til Solaris 2000 platforme

Til Solaris 4000 platforme

	Artikelnr.	Univ	ersal	Totrins	
		18 x 30	36 x 24	18 x 30	
Holder til firkantede flasker, 125 ml	75004106	28	54	56	
Holder til firkantede flasker, 250 ml	75004107	28	52	56	
Holder til firkantede flasker, 500 ml	75004108	15	24	30	
Holder til firkantede flasker, 1 000 ml	75004109	15	24	28	

Tabel 29: Tilgængelige holdere til firkantede beholdere til Solaris 4000 platforme

Til Solaris 2000 I / 2000 R platforme

	Artikelnr.	Universal
		11 x 14
Holder til firkantede flasker, 125 ml	75004106	9
Holder til firkantede flasker, 250 ml	75004107	7
Holder til firkantede flasker, 500 ml	75004108	3
Holder til firkantede flasker, 1 000 ml	75004109	2

Tabel 30: Tilgængelige holdere til firkantede beholdere til Solaris 2000 I / 2000 R platforme

Til Solaris 4000 I / 4000 R platforme

	Artikelnr.	Universal
		18 x 18
Holder til firkantede flasker, 125 ml	75004106	16
Holder til firkantede flasker, 250 ml	75004107	16
Holder til firkantede flasker, 500 ml	75004108	9
Holder til firkantede flasker, 1 000 ml	75004109	9

Tabel 31: Tilgængelige holdere til firkantede beholdere til Solaris 4000 I / 4000 R platforme

1. 2. 8. Nalgene bægerglasstativ

Til Solaris 2000 platforme



	Artikelnr.		Universal		Toti	rins
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Størrelse 11 x 14 / 12 x 14			St	ativer/pladse	er	
Nalgene 30 ml bægerglas, 36 pladser	75004116	1/36	-	-	2/72	-
Nalgene 50 ml bægerglas, 28 pladser	75004129	1 / 28	-	-	2/56	-
Nalgene 100 ml bægerglas, 16 pladser	75004118	1 / 16	-	-	2/32	-
Nalgene 250 ml bægerglas, 9 pladser	75004119	1/9	-	-	2/18	-
Nalgene 400 ml bægerglas, 8 pladser	75004120	1/8	-	-	2/16	-
Nalgene 600 ml bægerglas, 5 pladser	75004121	1/5	-	-	2/10	-
Størrelse 9 x 18			St	ativer/pladse	ər	
Nalgene 30 ml bægerglas, 32 pladser	75004110	-	2/64	2/64	-	4 / 128
Nalgene 50 ml bægerglas, 24 pladser	75004128	-	2/48	2 / 48	-	4 / 96
Nalgene 100 ml bægerglas, 15 pladser	75004112	-	2/30	2/30	-	4 / 60
Nalgene 250 ml bægerglas, 8 pladser	75004113	-	2/16	2/16	-	4/32
Nalgene 400 ml bægerglas, 6 pladser	75004114	-	2/12	2/12	-	4 / 24
Nalgene 600 ml bægerglas, 4 pladser	75004115	-	2/8	2/8	-	4 / 16

Tabel 32: Tilgængelige Nalgene bægerglasholdere til Solaris 2000 platforme

Til Solaris 4000 platforme

	Artikelnr.	Universal		Totrins	
		18 x 30	36 x 24	18 x 30	
Størrelse 11 x 14 / 12 x 14			Stativer/pladser		
Nalgene 30 ml bægerglas, 36 pladser	75004116	-	-	-	
Nalgene 50 ml bægerglas, 28 pladser	75004129	-	-	-	
Nalgene 100 ml bægerglas, 15 pladser	75004118	-	-	-	
Nalgene 250 ml bægerglas, 9 pladser	75004119	-	-	-	
Nalgene 400 ml bægerglas, 8 pladser	75004120	-	_	-	
Nalgene 600 ml bægerglas, 5 pladser	75004121	-	-	-	
Størrelse 9 x 18			Stativer/pladser		
Nalgene 30 ml bægerglas, 32 pladser	75004110	3 / 96	4 / 128	6 / 192	
Nalgene 50 ml bægerglas, 24 pladser	75004128	3/72	4 / 96	6/144	
Nalgene 100 ml bægerglas, 15 pladser	75004112	3 / 45	4 / 60	6 / 90	
Nalgene 250 ml bægerglas, 8 pladser	75004113	3/24	4 / 32	6 / 48	
Nalgene 400 ml bægerglas, 6 pladser	75004114	3 / 18	4 / 24	6/36	
Nalgene 600 ml bægerglas, 4 pladser	75004115	3/12	4 / 16	6/24	

Tabel 33: Tilgængelige Nalgene bægerglasholdere til Solaris 4000 platforme

Til Solaris 2000 I / 2000 R platforme

	Artikelnr.	Universal
Størrelse 11 x 14 / 12 x 14		Stativer/pladser
Nalgene 30 ml bægerglas, 36 pladser	75004116	1/36
Nalgene 50 ml bægerglas, 28 pladser	75004129	1 / 28
Nalgene 100 ml bægerglas, 16 pladser	75004118	1 / 16
Nalgene 250 ml bægerglas, 9 pladser	75004119	1/9
Nalgene 400 ml bægerglas, 8 pladser	75004120	1/8
Nalgene 600 ml bægerglas, 5 pladser	75004121	1/5
Størrelse 9 x 18		Stativer/pladser
Nalgene 30 ml bægerglas, 32 pladser	75004110	-
Nalgene 50 ml bægerglas, 24 pladser	75004128	-
Nalgene 100 ml bægerglas, 15 pladser	75004112	-
Nalgene 250 ml bægerglas, 8 pladser	75004113	-
Nalgene 400 ml bægerglas, 6 pladser	75004114	-
Nalgene 600 ml bægerglas, 4 pladser	75004115	-

Tabel 34: Tilgængelige Nalgene bægerglasstativer til Solaris 2000 I / 2000 R platforme
Til Solaris 4000 I / 4000 R platforme

	Artikelnr.	Universal
		18 x 18
Størrelse 11 x 14 / 12 x 14		Stativer/pladser
Nalgene 30 ml bægerglas, 36 pladser	75004116	-
Nalgene 50 ml bægerglas, 28 pladser	75004129	-
Nalgene 100 ml bægerglas, 16 pladser	75004118	-
Nalgene 250 ml bægerglas, 9 pladser	75004119	-
Nalgene 400 ml bægerglas, 8 pladser	75004120	-
Nalgene 600 ml bægerglas, 5 pladser	75004121	-
Størrelse 9 x 18		Stativer/pladser
Nalgene 30 ml bægerglas, 32 pladser	75004110	2 / 64
Nalgene 50 ml bægerglas, 24 pladser	75004128	2 / 48
Nalgene 100 ml bægerglas, 15 pladser	75004112	2 / 30
Nalgene 250 ml bægerglas, 8 pladser	75004113	2/16
Nalgene 400 ml bægerglas, 6 pladser	75004114	2/12
Nalgene 600 ml bægerglas, 4 pladser	75004115	2/8

Tabel 35: Tilgængelige Nalgene bægerglasstativer til Solaris 4000 I / 4000 R platforme

9

1. 2. 9. Skilletragtholder

Til Solaris 2000 platforme

	Artikelnr.	Universal			Toti	rins
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Enkelt kolbeholder til skilletragt, 250 ml til 2 000 ml	75004125	2	З	4	2	5
Skilletragtholder, monteret vertikalt		4	4	6	-	-

Tabel 36: Tilgængelige skilletragtholdere til Solaris 2000 platforme

Til Solaris 4000 platforme

	Artikelnr.	Universal		Totrins
		18 x 30	36 x 24	18 x 30
Enkelt kolbeholder til skilletragt, 250 ml til 2000 ml	75004125	5	7	9
Skilletragtholder, monteret vertikalt		8	12	-

Tabel 37: Tilgængelige skilletragtholdere til Solaris 4000 platforme

Til Solaris 2000 I / 2000 R platforme

	Artikelnr.	Universal
		11 x 14
Enkelt kolbeholder til skilletragt, 250 ml til 2000 ml	75004125	-
Skilletragtholder, monteret vertikalt		-

Tabel 38: Tilgængelige skilletragtholdere til Solaris 2000 I / 2000 R platforme

Til Solaris 4000 I / 4000 R platforme

	Artikelnr.	Universal
		18 x 18
Enkelt kolbeholder til skilletragt, 250 ml til 2 000 ml	75004125	-
Skilletragtholder, monteret vertikalt		-

 Tabel 39: Tilgængelige skilletragtholdere til Solaris 4000 I / 4000 R platforme

1. 2. 10. Universalbakke

Til Solaris 2000 platforme

Universalbakke

11 x 14 / 12 x 14 Universalbakke 9 x 18



2

З

4

Tabel 40: Tilgængelige universalbakker til Solaris 2000 platforme

75004122

Til Solaris 4000 platforme

	Artikelnr.	Universal		Totrins
		18 x 30	36 x 24	18 x 30
Universalbakke 11 x 14 / 12 x 14	75004123	-	-	-
Universalbakke 9 x 18	75004122	3	6	6

Tabel 41: Tilgængelige universalbakker til Solaris 4000 platforme

Til Solaris 2000 I / 2000 R platforme

	Artikelnr.	Universal
Universalbakke 11 x 14 / 12 x 14	75004123	1
Universalbakke 9 x 18	75004122	-

 Tabel 42: Tilgængelige universalbakker til Solaris 2000 I / 2000 R platforme

Til Solaris 4000 I / 4000 R platforme

	Artikelnr.	Universal
		18 x 18
Universalbakke 11 x 14 / 12 x 14	75004123	-
Universalbakke 9 x 18	75004122	2

 Tabel 43: Tilgængelige universalbakker til Solaris 4000 I / 4000 R platforme

Tilbehør til universalbakker

	Artikelnr.	Universal
Reserve-O-ringe (universalbakke)	75004132	-

Tabel 44: Tilgængeligt tilbehør til universalbakker

1. 2. 11. Klæbemåtter

Generelle oplysninger om valg af klæbemåtter:

Måtter med en størrelse på 170 x 280 mm egner sig optimalt til platforme på 11 x 14 og 12 x 14.



Alle måtter kan tilskæres til enhver ønsket størrelse.

Til Solaris 2000 platforme

	Artikelnr.	ι	Jniversa	I	Totr unive	ins, ersal
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Måtte med god vedhæftning, 200 x 200 mm, pakke: 2 stk.	75004126	1	4	6	2	8
Måtte med god vedhæftning, 280 x 170 mm, pakke: 2 stk.	75004127	2	2	4	4	4
Måtte med dårlig vedhæftning, 200 x 200 mm, pakke: 2 stk.	75004111	1	4	6	2	8
Måtte med dårlig vedhæftning, 280 x 170 mm, pakke: 2 stk.	75004117	2	2	4	4	4

Måtterne sælges i pakker á 2 stk. Antallet af måtter, der passer på en platform, relaterer ikke til denne pakkestørrelse, men til en enkelt måtte.

Tabel 45: Tilgængelige klæbemåtter til Solaris 2000 platforme

Til Solaris 4000 platforme

	Artikelnr.	Universal		Totrins	
		18 x 30	36 x 24	18 x 30	
Måtte med god vedhæftning, 200 x 200 mm, pakke: 2 stk.	75004126	6	12	12	
Måtte med god vedhæftning, 280 x 170 mm, pakke: 2 stk.	75004127	6	9	12	
Måtte med dårlig vedhæftning, 200 x 200 mm, pakke: 2 stk.	75004111	6	12	12	
Måtte med dårlig vedhæftning, 280 x 170 mm, pakke: 2 stk.	75004117	6	9	12	

Måtterne sælges i pakker á 2 stk. Antallet af måtter, der passer på en platform, relaterer ikke til denne pakkestørrelse, men til en enkelt måtte.

Tabel 46: Tilgængelige klæbemåtter til Solaris 4000 platforme

Til Solaris 2000 I / 2000 R platforme

	Artikelnr.	Universal
Måtte med god vedhæftning, 200 x 200 mm, pakke: 2 stk.	75004126	1
Måtte med god vedhæftning, 280 x 170 mm, pakke: 2 stk.	75004127	2
Måtte med dårlig vedhæftning, 200 x 200 mm, pakke: 2 stk.	75004111	1
Måtte med dårlig vedhæftning, 280 x 170 mm, pakke: 2 stk.	75004117	2

Måtterne sælges i pakker á 2 stk. Antallet af måtter, der passer på en platform, relaterer ikke til denne pakkestørrelse, men til en enkelt måtte.

Tabel 47: Tilgængelige klæbemåtter til Solaris 2000 I / 2000 R platforme

Til Solaris 4000 I / 4000 R platforme

	Artikelnr.	Universal
		18 x 18
Måtte med god vedhæftning, 200 x 200 mm, pakke: 2 stk.	75004126	4
Måtte med god vedhæftning, 280 x 170 mm, pakke: 2 stk.	75004127	2
Måtte med dårlig vedhæftning, 200 x 200 mm, pakke: 2 stk.	75004111	4
Måtte med dårlig vedhæftning, 280 x 170 mm, pakke: 2 stk.	75004117	2

Måtterne sælges i pakker á 2 stk. Antallet af måtter, der passer på en platform, relaterer ikke til denne pakkestørrelse, men til en enkelt måtte.

Tabel 48: Tilgængelige klæbemåtter til Solaris 4000 I / 4000 R platforme

1. 2. 12. Generelt tilbehør

1	. 2. 12. Generelt tilbehø	r	No.
		Artikelnr.	- V
	Applikatorrulle til klæbemåtter	75004124	
	Skruetrækker til tilbehørsdele	75004131	
	Gasfordeler (Solaris 2000 I / 2000 R)	SK2000-8GM	
	Gasfordeler (Solaris 4000 I / 4000 R)	SK4000-8GM	
	Kolbeholder-reservedelssæt (skruer)	SK0010	
	Platform-reservedelssæt til Solaris 2000, 2000 I/R og 4000 I/R (platformsskruer, værktøj)	SK0100	
	Platform-reservedelssæt til SK4000 (platformsskruer, værktøj)	SK0101	

Tabel 49: Tilgængeligt almindeligt tilbehør

1.3. Standarder og direktiver

Region	Direktiv	Standarder
Europa	2006/42/EF	EN 61010-1
	Maskindirektivet	IEC 61010-2-051
	2011/65/EF RoHS	EN 61326-1 Klasse B
	Begrænsning af anvendelsen af	EN ISO 14971
	bestemte farlige stoffer i elektrisk og elektronisk udstyr	EN ISO 9001
	Beskyttelsesmål:	Kun til Solaris 2000 I / 2000 R / 4000 I / 4000 B [,] IEC 61010-2-010
	2014/35/EF	<u>+00017 +00011.</u> IEO 01010 2 010
	Lavspændingsdirektivet	
	2014/30/EC Elektromagnetisk kompatibilitet (EMC)	
Nordamerika		ANSI/UL 61010-1
		IEC 61010-2-051
		IEC 61326-1 Klasse B
		CFR 47 FCC 15 EMC
		EN ISO 14971
		EN ISO 9001
		<u>Kun til Solaris 2000 I / 2000 R /</u> <u>4000 I / 4000 R:</u> IEC 61010-2-010
Japan		IEC 61010-1
		IEC 61010-2-051
		IEC 61326-1 Klasse B
		EN ISO 14971
		EN ISO 9001
		Kun til Solaris 2000 I / 2000 R / 4000 I / 4000 R: IEC 61010-2-010

 Tabel 50:
 Standarder og direktiver

BEMÆRK: Dette apparat er blevet testet og overholder grænseværdierne for et digitalt apparat klasse A Del 15 i FCC-bestemmelserne. Disse grænseværdier skal give en passende beskyttelse mod skadelig interferens, hvis apparatet anvendes i erhvervsmæssige omgivelser. Dette apparat genererer, anvender og kan afgive højfrekvent energi og kan, hvis det ikke installeres og anvendes i overensstemmelse med brugsanvisningen, forårsage skadelig interferens i radiokommunikation. Brugen af dette apparat i et boligområde kan medføre skadelig interferens. I dette tilfælde er brugeren forpligtet til at afhjælpe interferensen for egen regning.

1.4. Produktoversigt

1.4.1. Solaris 2000



① Platform-monteringsplade; ② Grafisk brugergrænseflade; ③ Hovedafbryder;

④ Lysnettilslutning; ⑤ Sikring; ⑥ Ethernet-tilslutning; ⑦ USB-tilslutning

Afbildning 1: Oversigt Solaris 2000

1.4.2. Solaris 4000



- ① Platform-monteringsplade; ② Grafisk brugergrænseflade; ③ Hovedafbryder;
- ④ Lysnettilslutning; ⑤ Sikring (kan nulstilles); ⑥ Ethernet-tilslutning; ⑦ USB-tilslutning

Afbildning 2: Oversigt Solaris 4000

Produktoversigt

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



① Platform-monteringsplade; ② Grafisk brugergrænseflade; ③ Hovedafbryder;

 \circledast Kabelgennemføring; \circledast Lysnettilslutning; \circledast Sikring (kan nulstilles); \oslash Ethernet-tilslutning;

⑧ USB-tilslutning; ⑨ Låg

Afbildning 3: Oversigt Solaris 2000 I / 2000 R

1.4.4. Solaris 4000 I / 4000 R



Afbildning 4: Oversigt Solaris 4000 I / 4000 R

1.4.5. Tilslutninger

Åbne rysteapparater



① Lysnettilslutning; ② Sikring; ③ Ethernet-tilslutning

Afbildning 5: Set bagfra (venstre: Solaris 4000, højre: Solaris 2000)

Inkuberede rysteapparater (kun opvarmede) og kølede rysteapparater (kølede og opvarmede)



1 Lysnettilslutning; 2 Sikring; 3 Ethernet-tilslutning; 4 Afløbsslange

Afbildning 6: Set bagfra (venstre: Solaris 2000 I / 2000 R, højre: Solaris 4000 I / 4000 R)

Lysnettilslutning

Rysteapparatet kræver en strømkilde, der svarer til dets specifikationer. Lysnetkabler medfølger.

Artikelnr.	Rysteapparat	Tekniske data
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

Afbildning 7: Specifikationer for rysteapparaters strømforsyning

ADVARSEL Beskadigelse på grund af forkert lysnettilslutning eller lysnetstik. Sørg for, at rysteapparatet kun tilsluttes stikkontakter, der er korrekt jordforbundet i overensstemmelse med de gældende elektriske sikkerhedsstandarder. Brug ikke rysteapparatet med et beskadiget eller utilstrækkelig dimensioneret lysnetkabel.

Netstikket skal til enhver tid være frit tilgængelig.

Sørg for, at området omkring rysteapparatet er tørt for at undgå risikoen for elektrisk stød. I tilfælde af utilsigtet spild af eller stænk fra væsker skal lysnetstikket trækkes ud af rysteapparatet, de spildte væsker skal fjernes, og enhver sundhedsmæssig eller biologisk fare skal elimineres, før du fortsætter.

Træk lysnetstikket til rysteapparatet ud, når det ikke er i brug.

<u>Sikring</u>

Sikringen kan nulstilles, hvis den er sprunget på grund af overstrøm. Rysteapparatet kan først tages i brug, når sikringen nulstilles i dens rigtige position.

Ethernet

Rysteapparatet er udstyret med en RJ45-Ethernet-tilslutning, der kan anvendes til at tilslutte det til et lokalt netværk (LAN). Der må kun anvendes apparater iht. standarden IEC 60950-1 med RJ45-Ethernet-tilslutning. Ved hjælp af J45-Ethernet-tilslutningen kan der downloades tilgængelige softwareopdateringer på et senere tidspunkt.

Kabelgennemføringer

Temperaturstyrede rysteapparater ha to åbninger i siden, der betegnes som kabelgennemføringer. Disse kabelgennemføringer er som regel lukket med kapper. Du kan fjerne kappen og føre ekstra apparater som f.eks. en ekstra temperatursensor ind eller føre kabler eller slanger som f.eks. gasfordelerens slanger igennem.

<u>USB</u>

Rysteapparatet har 2 USB-A 2.0-indgange, der kan anvendes til en almindelig USBnøgle. Der må kun anvendes apparater med USB-tilslutning iht. standarden IEC 60950-1.

Intern overtemperaturbeskyttelse

Temperaturstyrede rysteapparater har en termostat med manuel nulstilling, der er anbragt ved siden af varmeelementerne. Denne termiske anordning slukker for varmeelementerne for at undgå for høje temperaturer i rysteapparatets hus, hvis luftblandingsventilatoren skulle svigte.

Termostaten med manuel nulstilling skal nulstilles manuelt, for at rysteapparatet kan bruges igen. Dette kan kun gennemføres af Thermo Fisher Scientific-kundeservice.

2. Transport og opstilling



Kunden er selv ansvarlig for, at rysteapparatet opstilles korrekt.

BEMÆRK

Emballagen skal kontrolleres omgående ved levering. Kontroller emballagen omhyggeligt for transportskader før du pakker det leverede apparat ud. Hvis der konstateres en skade, skal fragtføreren notere skaden på din kopi af følgesedlen og underskrive den.

Åbn æsken forsigtigt og kontroller, at alle komponenter ("Tabel 51: Leveringsomfang") er leveret, før du bortskaffer emballagen. Hvis du konstaterer en skade efter udpakningen, skal du informere speditøren om dette og forlange en skadesundersøgelse.

Vigtigt: Hvis der ikke forlanges en skadesundersøgelse få dage efter modtagelsen af forsendelsen, fritages fragtføreren for erstatningsansvar. Du skal forlange en skadesundersøgelse.

2.1. Udpakning

Brug paklisten under udpakningen til at sikre at du har modtaget alle dele. Emballagen bør ikke bortskaffes, før det er kontrolleret, at alle dele er leveret.

Leveringsomfang

Artikel	Mængde
Rysteapparat	1
Lysnettilslutningskabel	1
Universalplatform	1*
Skruer til platform	
- Solaris 4000	4*
- andre versioner	3*
Brugsanvisning (print, en)	1
Brugsanvisning på USB-nøgle	1
Fastspændingsværktøj	1
(nøgle med T-greb)	
* På rysteapparaterne 2000 I, 2000 R, 40	00 I og 4000 R er platformen
allerede installeret i rysteapparatet. Hverke	en platformen eller skruerne fås
som enkeltdele.	

Tabel 51: Leveringsomfang

Kontakt Thermo Fisher Scientific, hvis du ikke har modtaget alle dele.

2.2. Opstillingssted

FORSIGTIG Beskyttelsen kan blive forringet af, at plastdele udsættes for ultraviolette stråler, og at deres stabilitet dermed reduceres. Udsæt ikke rysteapparatet og tilbehørsdele af plast for direkte sollys eller andre kilder af ultraviolette stråler.

Stil rysteapparatet på et plant bord eller en plan bænk, der kan bære rysteapparatets vægt med tilbehør og prøver, når det er i brug. Placer rysteapparatet i nærheden af en stikkontakt, der opfylder kravene på typeskiltet. Sørg for, at der er tilstrækkelig plads omkring apparatet til fri luftkonvektion, tilbehørsdele og nemmere betjening.

Følgende krav skal opfyldes ved opstilling af apparatet:

- For at sikre en tilstrækkelig afstand til bevægelige dele og undgå fastklemningsskader eller skader på apparater i nærheden skal der ved brug af et åbent rysteapparat være en afstand på 8 cm (3 in) på alle platformens fire sider. Det åbne rysteapparats platform kan rage ud over huset.
- Rysteapparater forårsager vibrationer. Der må ikke opbevares følsomme apparater eller farlige genstande eller stoffer i sikkerhedszonen.



Afbildning 8: Opstilling af åbne rysteapparater med en sikkerhedsafstand på 8 cm (3 in)

For at sikre en korrekt ventilation og en optimal ydelse skal følgende afstandskrav til temperaturstyrede rysteapparater overholdes:

- De inkuberede rysteapparater (kun opvarmede) Solaris 2000 I og 4000 I kræver en sikkerhedsafstand på 8 cm (3 in) på alle husets fire sider, se Afbildning 9.
- De kølede rysteapparater (kølede og opvarmede) Solaris 2000 R og 4000 R kan anvendes med en afstand på 8 cm (3 in) på alle husets fire sider, hvis der er et **frit område** som vist i venstre side i Afbildning 9.

 Hvis det frie område er monteret over det kølede rysteapparat (kølet og opvarmet) Solaris 2000 R og 4000 R, skal afstanden omkring huset forstørres til 30 cm (12 in) på alle fire sider som vist i højre side i Afbildning 9.



Afbildning 9: Opstillingsafstande for inkuberede rysteapparater (kun opvarmede) (venstre) og kølede rysteapparater (kølede og opvarmede) (højre) og kølede rysteapparater (kølede og opvarmede) med et frit område over apparater er vist i illustrationen til venstre.

ADVARSEL Under rysteprocessen skal farlige stoffer holdes borte fra denne zone. Underlaget skal:

- » være ren,
- » stabilt, robust, hårdt og vibrationsfrit,
- » give mulighed for at rysteapparatet kan opstilles horisontalt,
- » kunne bære rysteapparatets vægt.
- Rysteapparatet må kun anvendes indendørs.
- Rysteapparatet må hverken udsættes for varme eller kraftigt sollys.
- Opstillingsstedet skal altid være godt ventileret.
- Netstikket skal til enhver tid være frit tilgængelig.

2.3. Transport

FORSIGTIG Personskader på grund af at rysteapparatet falder ned. Løft altid rysteapparatet i begge sider. Løft aldrig rysteapparatet i frontpanelet eller i den installerede platform. Løft aldrig et temperaturstyret rysteapparat i det gennemsigtige lågs greb.

FORSIGTIG Rysteapparatets sikre drift påvirkes af kondensation af vand i huset. Dette sker, når rysteapparatet anbringes et varmere sted med en mere fugtig luft fra et køligere sted, f.eks. fra et lagerrum eller en transportvogn til laboratoriet. Lad apparatet tørre i mindst 2 timer før det tages i brug.

BEMÆRK Fjern altid åbne rysteapparatmodellers platform, deres fyldning og alle tilbehørsdele før du bevæger rysteapparatet. Hvis du ikke fjerner platformen, kan platform-monteringspladen eller rystemekanismen blive beskadiget. På temperaturstyrede rysteapparater er det nok af tage fyldningen af platformen.

BEMÆRK Apparatet kan blive beskadiget af stød. Transporter rysteapparatet i oprejst position og, om muligt, i den originale emballage.

Håndtering af rysteapparatet

Sørg under håndtering af rysteapparatet for:

- At løfte rysteapparatet i begge sider og aldrig foran og bag
- At der er så mange personer til at bære rysteapparatet, som dets vægt kræver, dog aldrig færre end to personer (du finder yderligere oplysninger i afsnittet "Tekniske specifikationer" på side 11).





Afbildning 10: Rysteapparatet bæres

ADVARSEL Løft altid rysteapparatet i begge sider. Løft aldrig rysteapparatet i for- eller bagsiden. Rysteapparaterne er tunge (du finder yderligere oplysninger i afsnittet "Tekniske specifikationer" på side 11). Det er nødvendigt, at mindst 2 personer løfter og bærer et rysteapparat.

FORSIGTIG Bær aldrig et temperaturstyret rysteapparat i det gennemsigtige lågs greb. Grebet kan brække af, og rysteapparatet kan falde på gulvet og kvæste personer alvorligt.

2.4. Nivellering

Selve rysteapparatet har ingen nivelleringsanordninger. Et nivelleret underlag er nødvendigt til korrekt opstilling af rysteapparatet.

BEMÆRK Læg ikke afstandsskiver eller andre flade genstande mellem rysteapparatets fødder og underlaget for at nivellere rysteapparatet.

2.5. Lysnettilslutning

Rysteapparatet kræver en strømkilde, der svarer til dets specifikationer. Lysnetkabler medfølger. Du finder yderligere oplysninger i Afbildning 7 på side 47.

ADVARSEL Beskadigelse på grund af forkert lysnettilslutning eller lysnetstik. Sørg for, at rysteapparatet kun tilsluttes til korrekt jordforbundne stikkontakter. Brug ikke rysteapparatet med et beskadiget eller utilstrækkelig dimensioneret lysnetkabel.

BEMÆRK Elektromagnetisk stråling kan medføre forstyrrelser på displayet. Dette beskadiger ikke apparatet og begrænser eller ændrer ikke dets funktion. For at undgå forstyrrelser på grund af elektromagnetisk stråling bør mobilenheder som f.eks. mobiltelefoner ikke anbringes i umiddelbar nærhed af apparatet. Brug ikke apparatet sammen med andre apparater med høj elektrisk effekt i en fælles strømkreds. Brug ikke flere apparater i en fælles multistikdåse. Sådan tilsluttes rysteapparatet til strømforsyningen:

- 1. Sluk apparatet med tænd-/sluk-kontakten på højre side.
- 2. Kontroller, at kablet overholder sikkerhedsreglerne i dit land.
- Sørg for, at netspænding og -frekvens stemmer overens med dataene på typeskiltet.

Netstikket skal til enhver tid være frit tilgængelig.

Sørg for, at området omkring rysteapparatet er tørt for at undgå risikoen for elektrisk stød. I tilfælde af utilsigtet spild af eller stænk fra væsker skal lysnetstikket til rysteapparatet trækkes ud, og de spildte væsker skal fjernes, før du fortsætter.

Træk lysnetstikket til rysteapparatet ud, når det ikke er i brug.

2.6. Grundindstilling

Før apparatet tages i brug, skal følgende trin gennemføres:

1. Tænd for apparatet. Thermo Scientific-logoet vises. Tryk på knappen Start Setup.



Afbildning 11: Opfordring til grundindstilling

2. Vælg det ønskede sprog i skærmbilledet Sprog. Tryk på knappen Næste.



Afbildning 12: Grundindstilling - Sprog

3. Du kan evt. indtaste apparatets navn i dialogboksen Apparatnavn. Tryk på knappen Næste.

	Unit N	ame	
Unit Name:	Tap to enter		
E	3ack	Next	

Afbildning 13: Grundindstilling - Apparatets betegnelse

 Du kan evt. indtaste by og land i tekstfeltet. Så snart du har indtastet tre tegn, bliver der i tekstfeltet By og land vist en liste med passende forslag. Du kan evt. vælge et forslag og derefter trykke på knappen Næste.

	Set Re	gion
	Enter the city and co	untry closest to your time zone
City, Country:	Tap to ente	r i

Afbildning 14: Grundindstilling - Region

5. Vælg det ønskede datoformat. Tryk på knappen Næste.

Set Date Di	isplay Format
(•) MN	MM/DD/YYYY
O DD,	/MMM/YYYY
Ow	YY/MMM/DD
Back	Next

Afbildning 15: Grundindstilling - Indstilling af datoformat

 Drej på hver drejekontakt på indstillingshjulet for at vælge den aktuelle dato. Tryk på knappen Næste.



Afbildning 16: Grundindstilling - Indstilling af dato

7. Vælg det ønskede klokkeslætformat og drej hver drejekontakt på indstillingshjulet for at indstille det aktuelle klokkeslæt. Tryk på knappen **Næste**.



Afbildning 17: Grundindstilling - Indstilling af klokkeslæt

 Kun inkuberede rysteapparater (kun opvarmede) og kølede rysteapparater (kølede og opvarmede): Valg af den ønskede temperaturenhed. Tryk på knappen Næste.

Temper	rature Units
	°C
C) 'F
Back	Next

Afbildning 18: Grundindstilling - Vælg temperaturenhed

 Kun inkuberede rysteapparater (kun opvarmede) og kølede rysteapparater (kølede og opvarmede): Klik på den blå pil for at indstille alarmen for overtemperatur og grænseværdierne for et nødstop. Tryk på knappen Næste.

High Temperature Alarm Settings	Low Temperature Alarm Settings				
High and Low Temperature Alarms will alert you if the shaker temperature deviates outside of the desired range. Default values are below, but you may change any of the settings.	High and Low Temperature Alarms will alert you if the shaker temperature deviates outside of the desired range. Default values are below, but you may change any of the settings.				
High Temp Alarm +5°℃ ➤	Low Temp Alarm -5°C 💙				
High Temp Shutdown +15°C 💙	Low Temp Shutdown -15°C 💙				
Back Next	Back Next				

Afbildning 19: Grundindstilling - Indstilling af alarmgrænseværdier for for høj og for lav temperatur

- Gentag fremgangsmåden for grænseværdierne for alarmen for undertemperatur. Tryk på knappen Næste for at fortsætte.
- Hvis du vil fortsætte uden at indstille en administrator-adgangskode, skal du lade valgmuligheden Uden adgangskontrol være valgt og trykke på knappen Næste for at fortsætte.

Hvis du vil indstille en administrator-adgangskode nu, skal du trykke på **Med** adgangskontrol.

Acces	ss Mode
Open Mode Select to allow full access	to all functionality for all users.
Secure Mode Select to require users to e changes to stored paramet (All users still able to use pro	nter the admin passcode to make ters, programs, and settings. ograms and set new run parameters,
Back	Next

Afbildning 20: Grundindstilling - Valg af adgangstilstand

12. Indtast en administrator-adgangskode i den viste opfordring til indtastning af en adgangskode med tastaturet og tryk på knappen Næste.



Afbildning 21: Grundindstilling - Valg af adgangstilstand

13. Indtast din administrator-adgangskode næste gang du bliver opfordret til at indtaste adgangskoden.

14. Indtast admin-adgangskoden igen som bekræftelse tredje gang du bliver opfordret til at indtaste adgangskoden.

	Adm	in Pass	scode			Adm	in Pass	code	
If you do not wi field	ish to change s below blank	the pa and ta	sscode at th p the Next b	his time, leave the putton.	F	Please re-ent	er the n	ew passcode.	
*	*	*	*	*	*	*	*	*	*
	7	8	9			7	8	9	
	4	5	6			4	5	6	
	1	2	3			1	2	3	
	0	Cle	ear			0	Cle	ear	

Afbildning 22: Grundindstilling - Indtastning og bekræftelse af en ny administratoradgangskode

15. Tryk på knappen **Næste** i det bekræftelsesskærmbillede, der vises, for at fortsætte.

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		

Afbildning 23: Grundindstilling - Bekræftelse af ændringen af administrator-

adgangskoden

16. Der vises et skærmbillede med generelle fysiske installationsanvisninger for rysteapparatet. Tryk på knappen Næste.

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		

Afbildning 24: Grundindstilling - Opstilling og tilslutning

17. Skærmbilledet Konfiguration afsluttet vises. Tryk på knappen Næste for at afslutte proceduren.

Setup Co	omplete
Setup is c	complete.
We recommend calibrat temperature for your appl calibrate the ten	ing the unit at the ideal ication. Would you like to nperature now?
Calibration can always be accesse	d under the main settings screen.
Calibrate	Finish

Afbildning 25: Grundindstilling - Kalibrering eller afslutning af grundindstillingen

BEMÆRK På inkuberede rysteapparater (kun opvarmede) og kølede rysteapparater (kølede og opvarmede) kan du trykke på **Kalibrer** og udføre proceduren, der beskrives i afsnittet "4. 6. Temperaturkalibrering" på side 127.

2.7. Opbevaring



Før rysteapparatet og tilbehørsdelene opbevares, skal hele systemet rengøres og, om nødvendigt, desinficeres eller dekontamineres. Opbevar ikke rysteapparatet og tilbehørsdelene

FORSIGTIG før kontamineringstilstanden entydigt er afklaret. Hvis du ikke er sikker på fremgangsmåden, bedes du kontakte Thermo Fisher Scientific-kundeservice ("Rengøring" på side 124, "Desinficering" på side 125 og "Dekontaminering" på side 125).

- Før rysteapparatet og tilbehørsdelene opbevares, skal de rengøres og efter behov desinficeres og dekontamineres.
- Før opbevaringen skal rysteapparatet og tilbehørsdelene være fuldstændig tørret.
- Opbevar rysteapparatet et tørt og støvfrit sted.
- Opbevar rysteapparatet stående på dets fødder.
- Opbevar ikke rysteapparatet i direkte sollys.

2.8. Forsendelse



FORSIGTIG

Før rysteapparatet og tilbehørsdelene sendes, skal hele systemet rengøres og, om nødvendigt, desinficeres eller dekontamineres. Opbevar ikke rysteapparatet og tilbehørsdelene før kontamineringstilstanden entydigt er afklaret. Hvis du ikke er sikker på fremgangsmåden, bedes du kontakte Thermo Fisher Scientific-kundeservice ("Rengøring" på side 124, "Desinficering" på side 125 og "Dekontaminering" på side 125).

Vær opmærksom på følgende før du sender rysteapparatet:

- Rysteapparatet skal være rengjort og dekontamineret.
- Dekontamineringen skal bekræftes med et dekontamineringscertifikat.
 Et dekontamineringscertifikat kan bestilles hos Thermo Fisher Scientifickundeservice.

3. Betjening

3.1. Tænd/sluk

Tryk på tænd-/sluk-kontakten på højre side for at tænde eller slukke rysteapparatet ((l) eller (0)).

Brugergrænsefladen viser Thermo Scientific logoet mens den booter.

Når brugergrænsefladen er klar, viser den rysteapparatets aktuelle driftstilstand.

3. 2. Grafisk brugergrænseflade

Startskærmbilledet er rysteapparatets grafiske brugergrænseflades standardskærmbillede. Det er dit udgangspunkt for brug af rysteapparatet.

På startskærmbilledet kan du:

- indstille de vigtigste driftsparametre som f.eks. rysteapparatets omdrejningstal, varighed og temperatur (kun temperaturstyrede rysteapparater)
- starte og stoppe rysteapparatet
- få vist oplysninger om driftstilstanden og redigere alarmmeldinger og advarsler
- navigere til andre skærmbilleder med oplysninger om driftstilstanden og indstillingsmulighederne

Skærmbilledets indhold i åbne rysteapparater og temperaturstyrede rysteapparater er en smule forskelligt.

I Afbildning 26 er der vist et eksempel på et åbent rysteapparats startskærmbillede.



Afbildning 26: Et åbent rysteapparats GUI-startskærmbillede

Et temperaturstyret rysteapparats startskærmbillede har ekstra temperaturreguleringer som vist i Afbildning 27.



Afbildning 27: Et temperaturstyret rysteapparats GUI-startskærmbillede

3. 2. 1. Indstilling af rysteapparatets vigtigste driftsparametre

Fra startskærmbilledet behøver du kun at trykke én gang for at komme til skærmbillederne med indstillingerne af rysteapparatets vigtigste driftsparametre.

Indstilling af omdrejningstal

1. Tryk på feltet **Faktisk omdrejningstal** (© i Afbildning 26 eller ⑦ i Afbildning 27) for at åbne skærmbilledet **Nominel værdi for omdrejningstal**, der er vist i Afbildning 28.



Afbildning 28: Indstillingsværdi for omdrejningstal

- Tryk på pilene over eller under hver drejekontakt på indstillingshjulet for at indstille det ønskede omdrejningstal. Alternativt kan du ændre det ønskede nominelle omdrejningstal ved at dreje på hver drejekontakt på indstillingshjulet enkeltvist.
- 3. Tryk på knappen **Gem**.

BEMÆRK Hvis et pop op-vindue vises, som informerer dig om, at den nominelle værdi for omdrejningstal er uden for det tilladte område, er din indstilling uden for det driftsomdrejningstalområde, der understøttes af rysteapparatet som angivet i afsnittet "Tekniske data" på side 11. Korriger din indstilling og fortsæt.

 Kvitter for den efterfølgende dialog Dine indstillinger blev gemt for at vende tilbage til hovedskærmbilledet.

Kalkulator til beregning af cirkulær bane: Beregning af omtrentligt omdrejningstal fra tidligere apparatindstilinger

I dialogboksen Nominel værdi for omdrejningstal kan du også trykke på **Beregn** omdrejningstal for at åbne skærmbilledet **Kalkulator til beregning af cirkulær bane**: Kalkulatoren til beregning af den cirkulære bane er nyttig, hvis du skifter fra et rysteapparat med en anden mekanisk opbygning til et rysteapparat i Solaris-serien, der er beskrevet i denne brugsanvisning. Denne kalkulator anvender forskellen i den cirkulære banes størrelse for at give dig et groft skøn over det nominelle omdrejningstal, der er nødvendigt for at opnå et lignende resultat.

BEMÆRK Anvend ikke resultatet, der er beregnet med kalkulatoren til beregning af den cirkulære bane til dine produktionsprøver med det samme, men gennemfør en række tests som bekræftelse. Cellerne kan alt efter de eksisterende forskydningskræfter vokse hurtigere eller langsommere, danne andre proteiner eller blive beskadiget (eller dø), fordi de fysiske betingelser ved skift til en anden cirkulær bane ændrer sig.

1. Vælg måleenheden (millimeter eller tommer).



Afbildning 29: Kalkulator til beregning af cirkulær bane

2. Vælg den forrige cirkulære banes størrelse. Tryk på knappen **Beregn**. Følgende vindue vises til beregning af omdrejningstallet.



Afbildning 30: Cirkulær bane beregnet

- 3. Tryk på knappen **OK**.
- Du kan beregne det foreslåede omdrejningstal for en kundespecifik cirkulær banes størrelse. For at udføre en beregning af en kundespecifik cirkulær bane skal du vælge Andre.
- 5. Indtast den cirkulære banes størrelse.



Afbildning 31: Beregning af en kundespecifik cirkulær banes størrelse

6. Tryk på knappen Beregn.

Vinduet Beregn omdrejningstal vises med omdrejningstallet for den indtastede størrelse.

Indstilling af driftstid

Rysteapparatet kan anvendes i kontinuerlig drift eller med en forudindstillet varighed. I kontinuerlig drift kan du efter ønske stoppe rysteapparatet manuelt. I drift med forudindstillet varighed stopper rysteapparatets drev automatisk, når timeren udløber. Sådan får du vist timeren:

- forløbet tid: hvor længe rysteapparatet har kørt, siden du har trykket på knappen Start, eller
- resterende tid: hvor længe rysteapparatet kører endnu, indtil timeren udløber

- 1. Tryk på feltet **Tid** (⑤ i Afbildning 26 og Afbildning 27) for at åbne skærmbilledet Tidstilstand, der er vist i Afbildning 32.
- 2. Vælg Drift med forudindstillet varighed.



Afbildning 32: Valg af tidstilstand

- 3. Indtast timerne og minutterne under Tidsindstilling.
- 4. Tryk på knappen **Gem**.
- For at vende tilbage til hovedskærmbilledet trykker du på den viste opfordring til kvittering **Dine indstillinger blev gemt**.

Indstilling af temperaturen

På temperaturstyrede rysteapparater kan du ved hjælp af feltet **Nominel temperaturværdi** til enhver tid indstille en temperatur til din anvendelse.

- 1. Tryk på feltet **Nominel temperaturværdi** (nederste område i pos. (6) i Afbildning 27), for at åbne skærmbilledet **Nominel temperaturværdi**, der er vist i Afbildning 33.
- Tryk på pilene over eller under hver drejekontakt på indstillingshjulet for at forudindstille standardtemperaturen som rysteapparatet skal bibeholde. Alternativt kan du ændre den ønskede temperatur ved at dreje på hver drejekontakt på indstillingshjulet enkeltvist.



Afbildning 33: Indstilling af den nominelle temperaturværdi: Skærmbillede med kalibreringsvisning

3. Tryk på knappen **Gem**.

BEMÆRK Hvis et pop op-vindue vises, som informerer dig om, at den nominelle temperaturværdi er uden for det tilladte område, er din indstilling uden for det driftstemperaturområde, der understøttes af rysteapparatet som angivet i afsnittet "Tekniske data" på side 11. Korriger din indstilling og fortsæt.

BEMÆRK Hvis et pop op-vindue vises, som informerer dig om, at den valgte temperatur er uden for det tilladte område på grund af den aktuelle omgivelsestemperatur, er omgivelsestemperaturen uden for rysteapparatets driftstemperaturområde som angivet i afsnittet "Tekniske data" på side 11. Tryk på knappen **OK** for at bekræfte at du har fået besked og for at fortsætte eller vælge en anden temperatur.

BEMÆRK Skærmbilledet **Nominel temperaturværdi** viser den kalibrerede temperatur og det offset, der er indstillet under kalibreringen som vist i afsnittet nederst til venstre i Afbildning 33. Kalibreringsprocessen beskrives i afsnittet "Temperaturkalibrering" på side 127.

 For at vende tilbage til hovedskærmbilledet trykker du på den viste opfordring til kvittering Dine indstillinger blev gemt.

BEMÆRK Temperaturstyrede rysteapparater reagerer på en bestemt måde, når den nominelle temperatur nås, hvilket betegnes som "oversvingning". En oversvingning af temperaturen betyder, at kammerets nominelle temperatur i første omgang overstyres til en lidt højere (eller lavere) temperatur og derefter nærmer sig den nominelle temperatur. Når den nominelle temperatur bliver tilnærmet, viser displayet ikke denne oversvingning. I stedet for vises kammertemperaturens stigning (eller fald), indtil den nominelle temperatur nås.

Sådan startes og stoppes rysteapparatet

1. Tryk på knappen **Start** for at starte rysteapparatet.

Knappen Start bliver til knappen Stop.



Afbildning 34: Knapperne Start hhv. Stop

2. Tryk på knappen Stop for at stoppe rysteapparatet.

3. 2. 2. Driftstilstand

Hvis rysteapparatet er i en god driftstilstand, vises der et grønt hjertesymbol på betjeningspanelet i området Driftstilstand og Statistiske data (7) i Afbildning 26 og (a) i Afbildning 27). Ved at trykke på det grønne hjertesymbol åbnes

skærmbilledet Driftstilstand. Skærmbilledet Driftstilstand indeholder statistiske data til brug af dit rysteapparat inkl. ryste- og driftstimer i alt og for sidste session. Temperaturstyrede rysteapparater viser desuden timerne for køle- og opvarmningsprocesser.

Status: Good		2	×	Status: Good		2	×
Total Run Time		Last Run Time Ses	ssion	Total Run Time		Last Run Time Ses	sion
Hours Shaking	232hrs	Hours Shaking	60brs	Hours Shaking	232hrs	Hours Shaking	60hrs
	2021110			Hours Powered	332hrs	Hours Powered	61hrs
Hours Powered	332hrs	Hours Powered	61hrs	Hrs tempering	168hrs	Hrs tempering	61hrs
			0				0

Afbildning 35: Rysteapparatets statistiske data: Åbent rysteapparat (venstre) og temperaturstyret rysteapparat (højre)

Advarsel

Når der udlæses en advarsel, viser betjeningspanelet en gul bjælke med rulletekst øverst i det aktuelle skærmbillede. Derudover aktiveres der en akustisk alarm. Den gule advarselsbjælke med rulletekst forsvinder, når meldingen er gået igennem to gange. Kun den gule advarselstrekant viser, at der

findes en eller flere advarsler for rysteapparatet. Advarselstrekanten har en blå cirkel med en hvid kant, hvor antallet af aktive advarsler vises. Ved at trykke på advarselstrekanten i området Info og Driftstilstand (Ø i Afbildning 26 og ® i Afbildning 27) åbnes der et skærmbillede med en liste over alle aktuelt aktive advarsler. Den nyeste advarsel vises udvidet, så du kan se alle oplysninger som vist i Afbildning 36. Du kan scrolle gennem listen og trykke på et vilkårligt punkt på listen for at udvide det og få vist flere informationer.



Afbildning 36: Advarselsliste

Du kan vælge den aktive advarsel ved at trykke på afkrydsningsfeltet ved siden af advarslen. Når du har trykket på knappen **Kvitter**, forsøger systemet at slette advarslen fra listen. Når alle advarsler er slettet, skifter symbolet Driftstilstand til det oprindelige grønne hjerte igen.

Alarmmelding

Hvis der udsendes en alarmmelding, stopper rysteapparatet omgående for at



forhindre skader på prøverne og/eller selve apparatet. Du skal kvittere for alarmmeldingen på betjeningspanelet, før du kan fortsætte med at bruge apparatet.

Når der udsendes en alarmmelding, viser betjeningspanelet en rød bjælke øverst i det aktuelle skærmbillede. I området Info & driftstilstand (⑦ i Afbildning 26 og (® i Afbildning 27) vises der en rød alarmklokke med lydbølger på begge sider. Derudover aktiveres der en konstant akustisk alarm.

Under den røde alarmbjælke viser en rullende tekstbesked en sammenfatning af den aktuelle alarmmelding. En Snooze-knap vises i højre side, med hvilken du midlertidigt kan gøre alarmmeldingen lydløs. Hvis årsagen til alarmen ikke afhjælpes i det tidsrum, alarmmeldingen er lydløs, aktiveres den akustiske alarm igen. Du kan vælge, hvor længe den skal være lydløs i indstillingerne. Yderligere oplysninger, se afsnittet "Alarmmeldinger/advarsler" på side 70.

Når du har trykket på klokkesymbolet i området Info & driftstilstand (⑦ i Afbildning 26 og ⑧ i Afbildning 27) vises der et skærmbillede, der viser fuldstændige detaljer om den aktuelt aktive alarmmelding, som vist i Afbildning 37. Du kan scrolle gennem listen og trykke på et vilkårligt punkt på listen for at udvide det og få vist flere informationer.



Afbildning 37: Alarmmeldingsliste

Du kan vælge den aktive alarmmelding ved at trykke på afkrydsningsfeltet ved siden af alarmen. Når du trykker på knappen **Kvitter**, forsøger systemet at slette alarmen fra listen, hvis årsagen er afhjulpet. Når alle alarmmeldinger er slettet, skifter symbolet Driftstilstand til det oprindelige grønne hjerte igen.

<u>Fejl</u>

Hvis der opstår en funktionsfejl, viser rysteapparatet en fejlmelding og stopper omgående for at undgå at beskadige prøverne og/eller selve apparatet. Skærmbilledet bliver rødt, og yderligere interaktion er ikke længere mulig. Der vises en fejlmelding sammen med en fejlkode som vist i eksemplet i Afbildning 38.



Afbildning 38: Skærmbillede med fejlmelding

Løs problemet på følgende måde og forsøg at genoptage driften:

- 1. Notér fejlkoden, der vises på skærmbilledet med fejlmeldingen, ned.
- 2. Tryk på knappen Kvitter for at gøre den akustiske alarm lydløs.
- 3. Start rysteapparatet igen ved at slukke for det og tænde det igen.
- Hvis fejlmeldingen fortsat vises, bedes du kontakte Thermo Fisher Scientifickundeservice og angive den fejlkode, der vises på skærmbilledet med fejlmeldingen.

3. 2. 3. Indstillinger

Det andet symbol i navigationslinjen er symbolet Indstillinger. Skærmbilledet nedenfor vises, når du trykker på **Indstillinger**. Skærmbilledet Indstillinger indeholder flere knapper, end der kan vises på én gang i displayet. Du skal rulle for at se resten af knapperne som vist i højre side i Afbildning 39.

	Shake Lab 1 11:22:24AM Settings		Shake Lab 1 11:22:24AM Settings	
¢.	Alarms and Alerts	î	Controls	î
	Access Control		Display	1
Ē	Calibration		📋 🔲 Files and Info	ł
	Controls	~	🐥 X Service 🖷	~
		0		0

Afbildning 39: Skærmbilledet Indstillinger

Alarmmeldinger/advarsler

På skærmbilledet Alarmmeldinger/advarsler kan du vælge, hvordan og hvornår alarmmeldinger og advarsler udlæses.



Afbildning 40: Valg af forudindstillinger for alarmmeldinger og advarsler

Lydstyrke: Med denne skyder kan du ændre lydstyrken for den lyd, der aktiveres ved en alarmmelding eller advarsel. Træk skyderen mod venstre for at reducere lydstyrken eller mod højre for at øge den og tryk derefter på **Gem**. Denne lydstyrkeindstilling vedrører både alarmmeldinger og advarsler.



Afbildning 41: Indstilling af lydstyrke for alarmmeldinger og advarsler

Alarm ved overtemperatur/alarm ved undertemperatur (kuntemperaturstyrede rysteapparater): Med disse to valgmuligheder kan du indstille en alarmgrænseværdi for over- og undertemperatur i forhold til den valgte nominelle temperaturværdi (se "Indstilling af temperaturen" på side 65). Så snart temperaturen i rysteapparatets kammer falder til under grænseværdien for undertemperatur eller stiger til over grænseværdien for overtemperatur, udsender rysteapparatet en alarmmelding. Vælg en temperaturgrænseværdi og tryk på **Gem**.



Afbildning 42: Indstilling af alarmgrænseværdier for for høj og for lav temperatur

BEMÆRK Alarmgrænseværdierne forskydes, når du ændrer den nominelle temperaturværdi.

Slukning ved overtemperatur/slukning ved undertemperatur (kun temperaturstyrede rysteapparater): Med disse to valgmuligheder kan du indstille en over- og undertemperaturgrænseværdi til sikkerhedsfrakobling i forhold til den valgte nominelle temperaturværdi (se "Indstilling af temperaturen" på side 65). Så snart temperaturen i rysteapparatets kammer falder til under grænseværdien for undertemperatur eller stiger til over grænseværdien for overtemperatur, slukker

rysteapparatet automatisk for at beskytte de behandlede prøver i kammeret. Vælg en grænseværdi og tryk på **Gem**.



Afbildning 43: Indstilling af over- og undertemperaturgrænseværdier for sikkerhedsfrakobling

BEMÆRK Fejlgrænseværdierne forskydes, når du ændrer den nominelle temperaturværdi.

Den lydløse tilstands varighed: Den lydløse tilstands varighed fastlægger, hvor længe en alarmmelding bliver gjort lydløs, når du trykker på knappen Snooze på den røde bjælke øverst i hovedvinduet (se "Alarmmelding" på side 68). Du kan indstille den lydløse tilstands varighed til 5 minutter, 10 minutter eller 15 minutter. Standardindstillingen er 10 minutter. Tryk på knappen **Gem** for at gemme ændringerne.



Afbildning 44: Indstilling af den lydløse tilstands varighed for alarmmeldinger

Deaktivering af kalibreringsmeldinger (kun temperaturstyrede rysteapparater): med disse valgmuligheder kan du deaktivere beskederne, der regelmæssigt påminder dig om at kalibrere temperaturmålingen på dit rysteapparat (du finder yderligere oplysninger i afsnittet "Temperaturkalibrering" på side 127). Du kan vælge, om du ønsker at deaktivere kalibreringspåmindelserne helt, eller kun mens et program udføres.

Disable Calibration Notifications Disable Calibration Notifications while a program is actively running

Afbildning 45: Deaktivering af kalibreringsmeldinger

Adgangskontrol

På skærmbilledet Adgangskontrol kan du tildele en administrator-adgangskode for at registrere processer med begrænsede adgangsrettigheder.

Rysteapparatet er som standard indstillet til drift uden adgangskontrol, dvs. at du ikke skal indtaste en adgangskode for at få adgang til apparatet og betjene det. Et skift til drift "Med adgangskontrol" forlanger en admin-adgangskode af hver bruger, der ønsker at ændre rysteapparatets indstillinger. Når valgmuligheden Med adgangskontrol er valgt, vises knappen Log på i hjørnet øverst til højre i touchscreendisplayet.

Sammenligningen af drift med eller uden adgangskontrol nedenfor viser, hvornår det er nødvendigt at indtaste en adgangskode.

Grafisk brugergrænseflade

Foranstaltning	Adgangskode nødvendig i tilstand uden adgangskontrol	Adgangskode nødvendig i tilstand med adgangskontrol
Indstilling af nominelle værdier for omdrejningstal, varighed og temperatur	Nej	Nej
Start og bremsning af rysteapparatet	Nej	Nej
Beregning af cirkulære baner	Nej	Nej
Kvittering for alarmmeldinger og advarsler	Nej	Nej
Visning af skærmbilledet Filer og info	Nej	Nej
Visning af driftstilstand	Nej	Nej
Visning og eksport af hændelseslog og diagrammer	Nej	Nej
Midlertidig deaktivering af alarmmeldingers lyd	Nej	Nej
Brug af programmer	Nej	Nej
Oprettelse, redigering eller sletning af programmer	Nej	Ja
Ændring af displayindstillinger	Nej	Ja
Ændring af konfigurationsindstillinger	Nej	Ja
Ændring af alarm- og advarselsindstillinger	Nej	Ja
Ændring af indstillingerne for adgangskontrol	Nej	Ja
Import og eksport af programmer	Nej	Ja
Nulstilling til fabriksindstilling	Nej	Ja
Gennemførelse af en kalibrering	Nej	Ja
Installation af firmwareopgradering	Ja	Ja

Tabel 52: Obligatorisk indtastning af en adgangskode ved drift med eller uden adgangskontrol
Grafisk brugergrænseflade

Afbildning 46 viser skærmbilledet Adgangskontrol uden og med adgangskontrol.

< *	Shake Lab 1 11:22:24AM Access Control		< *	Shake Lab 1 11:22:24AM Access Control	Logout Admin
¢.	 Open Secure Will require passcode a 	access to	۵. •	Open	le access to
	change certain parame	eters		change certain para	meters
,	Change Admin Passcode	>	,	Change Admin Passcode	>
	Save	0	0	Save	0

Afbildning 46: Adgangskontrol: Uden adgangskontrol (venstre) og Med adgangskontrol (højre)

Hvis du trykker på **Med adgangskontrol** og derefter på **Gem**, bliver du opfordret til at indtaste admin-adgangskoden, før ændringen kan bekræftes. Alle apparater leveres med den samme adgangskode for administratorer, der er forprogrammeret på fabrikken. Adgangskoden er trykt i håndbogen.

Enhver bruger, der forsøger at ændre en indstilling, der er beskyttet med en adgangskode, bliver også opfordret til at indtaste admin-adgangskoden.

BEMÆRK Adgangskoden 00000 er fabriksindstillet for administratoren.

Med knappen **Skift admin-adgangskode** kan du ændre den forudindstillede adgangskode. Hvis du trykker på knappen **Skift admin-adgangskode**, bliver du opfordret til først at indtaste den aktuelle adgangskode og derefter den nye adgangskode og til sidst den nye adgangskode igen som bekræftelse. Beskeden **Din nye adgangskode blev gemt** viser, at ændringen blev udført uden fejl.

Kalibrering

Skærmbilledet Kalibrering (kun temperaturstyrede rysteapparater) åbner en række skærmbilleder, med hvilke du kan kalibrere rysteapparatet til en præcis temperaturstyring. Denne procedure beskrives i afsnittet "Temperaturkalibrering" på side 127.

Driftsparametre

Skærmbilledet Driftsparametre har valgmuligheder til forudindstilling af standarddriftsparametrene, med hvilke apparatet skal starte:

Indstillingsværdi for omdrejningstal: Du kan forudindstille den nominelle værdi for omdrejningstal til en værdi mellem 15 og 525 o/min. Drej på drejekontakterne på indstillingshjulet for at indstille omdrejningstallet og tryk på knappen **Gem**. Du finder detaljerede anvisninger om anvendelse af skærmbilledet Nominel værdi for omdrejningstal i afsnittet "Indstilling af omdrejningstal" på side 62.

Tidstilstand: Rysteapparatet kan anvendes i kontinuerlig drift eller med en forudindstillet varighed. I kontinuerlig drift kan du efter ønske stoppe rysteapparatet manuelt. I drift med forudindstillet varighed stopper rysteapparatets drev automatisk, når timeren udløber. Til drift med forudindstillet varighed kan du specificere en standarddriftstid og vælge, om brugergrænsefladen skal vise den forløbne eller den resterende tid, mens rysteapparatet kører. Du finder detaljerede anvisninger om anvendelse af skærmbilledet Tidstilstand i afsnittet "Indstilling af driftstid" på side 64.

Nominel temperaturværdi (kun temperaturstyrede rysteapparater): Med denne valgmulighed kan du indstille en standardtemperatur, som rysteapparatet skal vise, når det starter. Du finder detaljerede anvisninger om anvendelse af skærmbilledet Nominel temperaturværdi i afsnittet "Indstilling af temperaturen" på side 65.

Automatisk genstart: Med denne funktion genstartes apparatet efter et strømsvigt, der er opstået under et normalt start/stop-forløb, udførelse af et program eller en temperaturkalibrering. Hvis valgmuligheden Automatisk genstart er indstillet til **Nej**, genstarter apparatet ikke ved et strømsvigt.

<u>Display</u>

I Display-indstillingerne kan du ændre forskellige display-valgmuligheder.

Lysstyrke: Brug skyderen eller knapperne + / - til at indstille displayets lysstyrke.



Afbildning 47: Indstilling af skærmens lysstyrke

Sprog: Tryk på knappen **Sprog** for at ændre displayets sprog. Vælg det ønskede sprog ved at dreje på indstillingshjulet og tryk på knappen **Gem**.



Afbildning 48: Valg af displaysprog

Måleenheder (kun temperaturstyrede rysteapparater): Tryk på Enheder for at skifte alle temperaturvisninger på GUI-skærmbilleder mellem Grader celsius °C og Fahrenheit °F.

Units of Measure	°C 🔨
	°F

Afbildning 49: Valg af temperaturvisningsenhed

Dato: Tryk på knappen Dato for at indstille datoen. Drej på drejekontakterne på indstillingshjulet for måned, dag og år og tryk på knappen **Gem**.



Afbildning 50: Indstilling af måned, dato og klokkeslæt

Datoformat: Tryk på knappen **Datoformat** for at indstille datoformatet. Tryk på afkrydsningsfeltet for det ønskede datoformat (eksempel: MM/DD/ÅÅÅÅ vises som april 15, 2015) og tryk på knappen **Gem**.



Afbildning 51: Indstilling af datoformat

Tid: Tryk på knappen **Klokkeslæt** for at indstille tiden og tidsformatet. Tryk på **12 timer** eller **24 timer**, drej på drejekontakterne på indstillingshjulet for timer, minutter og AM/FM (kun 12-timers format) og tryk på knappen **Gem**.





Dvaletilstand: Tryk på **Dvaletilstand** så rysteapparatets display skifter til dvaletilstand efter 15 minutters inaktivitet. I dvaletilstand viser skærmbilledet opfordringen **Berør** for at aktivere.



Afbildning 53: Aktivering af dvaletilstand

Region: Tryk på knappen **Region** for at indstille den region, hvor apparatet anvendes. Tryk på tekstfeltet By, Land og begynd med at indtaste de første bogstaver i din bys navn. Brugergrænsefladen foreslår mulige overensstemmelser efter de første tre tegn. Accepter en overensstemmelse eller fortsæt med at indtaste det fulde navn og tryk derefter på **Gem** på displaytastaturet.

< *	Shi 11: Rei	ake La 22:24) gion	b 1 AM	Enter the	city and		closest to	your tim	ie zone		Shake Lab 1 11:22:24AM Region	
≎ ₀ <u>≅</u>	Cit	y, Coui	ntry:	N							Enter the city and country closest to your time City, Country: New York, United States	e zone
Q	w	E	R	Ţ	Y	U	Ţ	0	Ρ			
A	S	D	F	G	H	J	K	L	S	ave		
1	z	x	С	V	В	N	м			1		
?!1:	23								-		Save	0

Afbildning 54: Indstilling af region

Apparatbetegnelse: Tryk på knappen Apparatnavn for at navngive eller omdøbe rysteapparatet. Tryk på tekstfeltet Apparatnavn og begynd med indtastningen. Tryk på Gem på displaytastaturet når du er færdig.

< ♠	Sh 11 Ur	ake La :22:24) hit Nam	b 1 AM ie			2					< ☆	Shake Lab 1 11:22:24AM Unit Name	Ø	
٥.	U	nit Na	me	Sha	ake La	ab 1]	۵.			
Q	W	E	R	т	Y	U	T	0	Р			Unit Name:	Shake Lab 1	
А	S	D	F	G	Н	J	К	L	Sa	ive				
1	Z	X	С	۷	В	Ν	М	@	+	1	÷			
?!1	23							-	-	.			Save	0

Afbildning 55: Navngivelse af rysteapparatet

Menuen Kundespecifik tilpasning: Tryk på knappen Menuen Kundespecifik tilpasning for at tilpasse de nederste to symboler på hoved-navigationslinjen kundespecifikt. Træk det ønskede symbol ud af hovedskærmbilledets område til det symbol du vil udskifte. Tryk på knappen **Gem** for at bekræfte.

BEMÆRK Ved at trykke på **Menuen Nulstilling** kan du til enhver tid nulstille navigationslinjen til fabriksindstillinger.



Afbildning 56: Tilpasning af navigationslinjen

Filer og info

Dette skærmbillede viser serienummeret, GUI-programsoftwaren og dens version, rysteapparatets hovedstyringsfirmware og dens version, parameterfilen og dens version og det aktuelle styresystem.

Du kan også trykke på **Nulstil til fabriksindstillinger** for at nulstille rysteapparatet til fabriksindstillingerne. Nulstilling til fabriksindstillingerne kræver en admin-adgangskode og sletter alle indstillinger, der er foretaget via den grafiske brugergrænseflade. Hændelsesprotokollen bliver ikke slettet ved nulstilling til fabriksindstillingerne.

<	Shake Lab 1 11.22:24AM Files and Info Serial Number 123TSX990960 H.M.N. 4.1.0 Main Controller 7000 0612 V3 Parameters 50154439 V1	î	<	Shake Lab 1 11:22:24AM Files and Info Main Controller 7000 0612 V3 Parameters 50154439 V1 Board Support Package: 4.1.15 jethro-15.0		^
Ē	Board Support Package: 4.1.15-jethro-15.0		Ē	Factory Reset	>	1
Ļ	Factory Reset	> 🗸	÷	Firmware	>	Ŷ
		()				()

Afbildning 57: Filer og info

Hvis du ruller længere ned, vises knappen **Firmware** som vist i højre side i Afbildning 57. Du kan trykke på denne knap for at installere en ny firmware til rysteapparatet. Du finder detaljerede anvisninger om installation af en ny firmware i afsnittet "4. 7. Firmwareinstallation" på side 131.

BEMÆRK En ny firmware skal købes hos en autoriseret servicetekniker.

<u>Service</u>

Adgangen til serviceindstillingerne er begrænset og er kun tilladt for autoriserede serviceteknikere.

3. 2. 4. Programmer

Tryk på symbolet **Programmer** på navigationslinjen for at få vist listen over programmer eller oprette et nyt program. Du kan oprette, redigere, slette, importere og eksportere et program. I Afbildning 58 er skærmbilledet Programmer vist for et nyt apparat (venstre) og et apparat, der har været anvendt i et stykke tid (højre). Det anvendte apparat viser en liste over de programmer, som brugeren har oprettet.



Afbildning 58: Programmer

Hvert program på programlisten har et øjesymbol **()**, der åbner en hurtig programvisning med detaljerede oplysninger om programmet og et blyantsymbol **/**, der åbner programmet til redigering.

Valg af forudindstillinger for alle programmer

Fra skærmbilledet Programmer kommer du ved hjælp af knappen **Udv. indstillinger** til skærmbilledet **Udvidede indstillinger**. På dette skærmbillede kan du vælge, hvad der sker, når et program startes. Disse indstillinger gælder for alle programmer, du opretter.



Afbildning 59: Indstilling af programspecifikationer

- 1. Tryk på knappen Udv. indstillinger.
- 2. Vælg en valgmulighed for **Vis tidsstyring** for at vælge, hvordan programtidsstyringen skal vises:
 - » Udløbet: hvor længe rysteapparatet har kørt, siden du har trykket på knappen Start, eller
 - » Resterende driftstid: hvor længe rysteapparatet kører endnu, indtil timeren udløber
- Hvis du ønsker, at rysteapparatet omgående starter, når et program startes, kan du aktivere valgmuligheden Start rysteapparat ved indlæsning af programmet. Som standard skal du trykke på knappen Start for at starte et program.

BEMÆRK Valgmuligheden **Start rysteapparatet ved indlæsning af programmet** fungerer kun som forventet, hvis rysteapparatets låg er lukket. Forløbet starter automatisk, så snart låget er lukket.

4. Tryk på knappen Gem.

Oprettelse af et program

Du kan oprette og gemme op til 99 programmer.

- 1. Tryk på knappen Opret nyt program.
- 2. Indtast programmets navn.



Afbildning 60: Skærmbilledet Oprettelse af program

3. Indtast programmets temperatur (kun temperaturstyrede rysteapparater), omdrejningstal og driftstid i timer og minutter.

Lad feltet **Temp** være tomt for at deaktivere temperaturstyringen på et temperaturstyret rysteapparat.

- For at tilføje et ekstra trin i dit program skal du rulle nedad og trykke på knappen Tilføj trin.
- 5. Tryk på knappen **Gem** for at gemme programmet.

Redigering af et program

- 1. Tryk på blyantsymbolet 🖍 ved siden af det program du ønsker at redigere.
- Rediger de ønskede felter. Tryk på knappen Gem. Programmet gemmes med de nye ændringer.
- Du kan tilføje et trin ved at vælge Tilføj trin nederst i skærmbilledet. Rul skærmbilledet nedad for at kunne se knappen Tilføj trin, hvis der findes flere end 3 trin.

BEMÆRK Hvis der vises en alarmmelding eller en fejlmelding under rysteprocessen med flere trin, stopper rysteapparatet automatisk. Hvis der vises en advarsel, fortsættes rysteprocessen.



Afbildning 61: Tilføj trin

Sletning af et program

- 1. Tryk på blyantsymbolet ved siden af det program du ønsker at redigere.
- 2. Tryk på knappen **Slet**. Der vises et vindue, hvor du anmodes om en bekræftelse.



Afbildning 62: Slet program

3. Tryk på knappen Slet. Det valgte program slettes.

Udførelse af et program

- 1. For at udføre et eksisterende program skal du vælge det program, du ønsker at udføre.
- Hvis du ønsker at se den hurtige visning af programmet, skal du trykke på øjesymbolet
 ved siden af programmet.
- 3. Vinduet Hurtig visning af program vises.



Afbildning 63: Hurtig visning af program

4. Tryk på **Run** for at indlæse programmet i hovedskærmbilledet.

Afhængigt af dine valgte **Automatisk drift**-indstillinger starter programmet omgående efter, du har trykket på knappen **Start** på startskærmbilledet.

 Når programmet er fuldstændig udført, vises følgende vindue. Tryk på knappen OK.



Afbildning 64: Program afsluttet

Import af programmer

Du kan importere programmer på et rysteapparat, der er oprettet på et andet. For at gøre dette skal USB-nøglen være sat i rysteapparatets USB-indgang.

Følgende skærmbillede viser alle programmer, der er identificeret på USB-nøglen.



Afbildning 65: Vælg de programmer, der skal importeres

Vælg de programmer du ønsker at importere. Tryk på knappen Import.

Når programmerne er blevet importeret, vises følgende vindue. Tryk på knappen **OK**. USB-nøglen kan nu fjernes.



Afbildning 66: Import afsluttet

Eksport af programmer

Du kan eksportere programmer fra et rysteapparat til et andet. Kontroller, at der er sat en USB-nøgle i.

Hvis USB-nøglen er sat i, vises følgende skærmbillede:



Afbildning 67: Vælg de programmer, der skal eksporteres

Vælg de programmer du ønsker at eksportere. Tryk på knappen Eksport.

Når eksporten er afsluttet, vises følgende vindue. Tryk på knappen **OK**. USB-nøglen kan nu fjernes.



Afbildning 68: Eksport afsluttet

<u>Hændelseslog</u>

Det tredje symbol i navigationslinjen er hændelsesloggen, som indeholder et datasæt med bruger- og systemhændelser. Skærmbilledet Hændelsesprotokol vises, så snart du trykker på symbolet for hændelseslog på navigationslinjen som vist i Afbildning 69.



Afbildning 69: Hændelseslog

På dette skærmbillede vises der en liste over alle hændelser, herunder dato og klokkeslæt.

Der kan vises yderligere oplysninger om en enkelt hændelse ved at vælge elementet på listen.

Hændelserne kan filtreres efter følgende kategorier:

- Alarmmeldinger
- Advarsler
- Ændringer af indstillingerne
- Start og stop
- Programkørsler
- Åbninger af låget (kun temperaturstyrede rysteapparater)

Hvis et filter er valgt, skifter knappen Vis alle til højre til Filter AKTIVERET som vist i Afbildning 70.

Event Log			
Date/Time	Event	Filter ON	~
5/12/15 11:34:06AM	Setpoint Change		~
5/12/15 11:34:06AM	High Temp Alarm		~
5/12/15 11:34:06AM	Unit Stopped		~
5/12/15	Unit Started		~

Afbildning 70: Filtreret hændelseslog

Eksport af hændelsesprotokol

1. Vælg de hændelser, der skal eksporteres, fra drop down-menuen **Hændelser, der skal eksporteres**. Vælg loggens eller rapportens eksportformat.



Afbildning 71: Valg af hændelseslog, der skal eksporteres

2. Der kan vælges et på forhånd defineret eller kundespecifikt datoområde.

BEMÆRK Du kan kun eksportere hændelser fra de sidste seks måneder.

 Eksporten kan gennemføres ved hjælp af en USB-nøgle. For at gemme loggen eller rapporten skal en USB-nøgle være sat i. Tryk på knappen Eksporter log for at downloade loggen eller rapporten.



Afbildning 72: Sæt en USB-nøgle i til eksporten

Diagrammer

Diagrammer viser data for omdrejningstal eller temperatur (kun temperaturstyrede rysteapparater) ved hjælp af tiden. X-aksen viser tiden og Y-aksen omdrejningstallet eller temperaturen.

Med temperaturdiagrammet kan du registrere den nominelle, faktiske temperatur og omgivelsestemperaturen tidsmæssigt. Med knappen **Rediger** kan du vælge, hvilke af de tre kurver, der skal optages i temperaturdiagrammet. Derudover kan visningsområdet ændres til 1 dag, 7 dage eller en brugerdefineret rækkefølge af dage, enten ved hjælp af alle 24 timer eller i valgte timer hver dag.

Knappen **Opdater** opdaterer diagramvisningen i overensstemmelse med dine indstillinger og viser alle nye data, der er registreret siden sidste indlæsning eller opdatering af diagrammet.

Showing	5/12/15, BAM 6	4PM	-	Se Se	tpoint	Actual		Showing: 5/	2/15, 8AM	to 4PM	Set	point	Actual	Ambie
350rpm							-	70.0°C						2
300rpm							¥0	60.0°C						
250rpm								50.0°C						
200rpm								40.0°C						
150rpm	-							30.0°C	-				1	
100rpm								20.0°C						
50rpm							1000	10.0°C						
Orom								0.0°C		0.114	10414	11444	100	

Afbildning 73: Omdrejningstal- og temperaturdiagrammer

Du kan sætte en USB-nøgle i rysteapparatets USB-indgang og trykke på **Eksporter diagramdata** for at downloade dataene for omdrejningstal- eller temperaturdiagrammet.

Redigering af diagrammet

1. Tryk på knappen **Rediger** for at begynde med redigeringen af diagrammet.



Afbildning 74: Redigering af diagrammet: Valg af temperatur eller omdrejningstal til registreringen

- Tryk på knappen Temp eller Omdrejningstal for at vælge, hvilket diagram, der skal vises.
- 3. Rul længere ned for at få vist valgmulighederne **Datoområde**.
- 4. Tryk på 1 dag, 7 dage eller **Brugerdefineret** for at vælge et brugerdefineret datoområde.
- 5. Drej på drejekontakterne på indstillingshjulet for at vælge startdato og tryk på knappen **Fra**.

 Drej på drejekontakterne på indstillingshjulet for at vælge slutdato og tryk på knappen Til.



Afbildning 75: Valg af datoområde til diagrammet

- 7. Rul længere ned for at få vist valgmulighederne Tidsperiode.
- 8. Tryk på 24 timer dagligt eller på knappen **Brugerdefineret** for at vise en bestemt periode og
 - a. Drej på drejekontakterne på indstillingshjulet for at vælge det daglige starttidspunkt og tryk på knappen **Fra**.
 - b. Drej på drejekontakterne på indstillingshjulet for at vælge det daglige sluttidspunkt og tryk på knappen **Til**.

BEMÆRK Ved at vælge start- og sluttidspunkterne vælger du en nøjagtig visning af perioden mellem to tidspunkter. Hvis du arbejder med valgmuligheden 24 timer dagligt, begynder registreringen kl. 00:00 og slutter kl. 23:59 hver dag.



Afbildning 76: Valg af tidspunkter på dagen til registreringen

9. Tryk på knappen **Gem** for at gemme ændringerne eller kundespecifikke indstillinger.

3.3. Tilbehør



Personskader og biologisk risiko på grund af ødelagte beholdere.

FORSIGTIG

Forkert installerede tilbehørsdele kan medføre glas, der går i stykker, og spildte prøver.

Kontroller, at tilbehørsdelene er installeret korrekt ved at anvende de rigtige værktøjer og skruer.

Kontroller, at tilbehørsdelene er placeret korrekt på platformen. Anvend altid kun beholdere med tilbehør, der er tilpasset den pågældende størrelse.



Risiko for snitsår på grund af skarpe kanter.

Vær forsigtig når du håndterer platforme og andre tilbehørsdele.

FORSIGTIG



Fingrene kan komme i klemme i en platform, der bevæger sig.

FORSIGTIG Læg aldrig fingrene på eller under en platform, der bevæger sig.

BEMÆRK

Kunden er selv ansvarlig for, at tilbehørsdelene installeres korrekt.

Den fremgangsmåde til installationen, der er beskrevet her, er indtil videre gældende for alle rysteapparat-udførelser, der er angivet i denne brugsanvisning.

Anvend altid de rigtige værktøjer, især de værktøjer, der er leveret med tilbehøret. Hvis du mister nogle af disse artikler, kan du efterbestille de reservedelssæt, der er angivet på listen ("Tilbehør" på side 17). Anvend kun det angivne værktøj og kun de medfølgende skruer.

3. 3. 1. Montering af platformen

FORSIGTIG Du finder en fuldstændig liste over kompatible platforme til hvert rysteapparat i afsnittet "1. 2. 1. Platforme" på side 18. Temperaturstyrede rysteapparater må kun anvendes med de platforme, der er forudinstalleret på fabrikken.

FORSIGTIG Risiko for snitsår på grund af skarpe kanter. Når du afmonterer en platform, skal du tage fat under platformen. Løft aldrig platformen i kolbeholderne, der er fastgjort til den. Vær forsigtig når du håndterer platforme og andre tilbehørsdele.

FORSIGTIG I rysteapparatet Solaris 4000 I / 4000 R er der en risiko for at få fingre og hænder i klemme, når 11 x 14-platformen bevæger sig. Anvend ikke 11 x 14-platformen til rysteapparatet Solaris 4000 I / 4000 R.

FORSIGTIG Anvend udelukkende det skruemateriale og fastspændingsværktøj, der medfølger til platformen. Anvendelse af andre skruer eller forkerte fastspændingsværktøjer medfører en forkert installation og dermed muligvis en beskadigelse af rysteapparatet og tilbehørsdelene.

Monter altid en platform med samtlige skruer.

Anvend altid det rigtige fastspændingsværktøj:

- 3/16"-fastspændingsværktøj (GT530066) til alle plafforme til rysteapparatet Solaris 2000, Solaris 2000 I/R og 4000 I/R
- 7/32"-fastspændingsværktøj (GT530080) til alle platforme til rysteapparatet Solaris 4000

<u>Universalplatforme</u>

Hvert rysteapparat i Solaris 2000- og 4000-serien leveres med en universalplatform, skruer og et fastspændingsværktøj. Du kan købe ekstra platforme til dine anvendelser. Du finder en fuldstændig liste over kompatible platforme til hvert rysteapparat i afsnittet "1. 2. 1. Platforme" på side 18.

På rysteapparaterne 2000 I, 2000 R, 4000 I og 4000 R er platformen allerede installeret i rysteapparatet. Hverken platformen eller skruerne fås som enkeltdele.



Afbildning 78: Universalplatform monteret på et Solaris 2000 rysteapparatet (venstre) og på et Solaris 4000 rysteapparat (højre)

 Læg platformen forsigtigt vandret på rysteapparatet og dets 3 monteringspunkter. Platforme til rysteapparatet Solaris 2000 har 3 monteringshuller. Platforme til rysteapparatet Solaris 4000 har 4 monteringshuller.

Platforme til rysteapparatet Solaris 2000 I/R og 4000 I/R har 3 monteringshuller.

 Spænd skruerne forsigtigt for at fastgøre platformen på rysteapparatet. Hvis fastspændingsværktøjet begynder at bøje, skal du stoppe med at spænde skruen.

Tilbehør

Totrins platforme

BEMÆRK Totrins platforme er ikke egnet til rysteapparatet Solaris 2000 I/R og 4000 I/R.



Afbildning 79: Totrins platformenhed

Forbind den øverste og nederste platform ved at montere dem med de 4 støtter i hvert hjørne. Monter støtterne med de pågældende skruer fra den øverste og den nederste platform.

Spænd skruerne forsigtigt for at forbinde støtterne og platformene. Hvis fastspændingsværktøjet begynder at bøje, skal du stoppe med at spænde skruen.

Den nederste platform indeholder monteringshullerne til fastgørelse på rysteapparatet. Kontroller, at den nederste platform befinder sig på den nederste side under samlingen.

Tilbehør

BEMÆRK Før du stiller beholdere på platformenheden, skal du gennemføre en sidste kontrol for at sikre, at platformenheden ikke vakler.



Afbildning 80: Totrins platform på et Solaris 4000 rysteapparat

1. Læg den monterede totrins platform forsigtigt vandret på rysteapparatet og dets platform-monteringsplade.

Platforme til Solaris 2000 har 3 monteringshuller, platforme til Solaris 4000 har 4.

 Spænd skruerne forsigtigt for at fastgøre platformen på rysteapparatet. Hvis fastspændingsværktøjet begynder at bøje, skal du stoppe med at spænde skruen.

3. 3. 2. Isætning af kolbeholder og beholder

FORSIGTIG Biologisk risiko på grund af ødelagte eller utætte beholdere. Forkert installerede tilbehørsdele kan medføre spildte prøver. Kontroller, at tilbehørsdelene er installeret korrekt ved at anvende de rigtige værktøjer og skruer. Til isætningen er en skruetrækker (PH2) med en skaftlængde på 150 mm/6 in nødvendig. (Artikelnummer 75004131). Kontroller, at tilbehørsdelene er placeret korrekt på platformen. Anvend altid en beholder med tilbehør, der er tilpasset den pågældende størrelse. Beholderne skal være intakte og installeret korrekt.

FORSIGTIG Risiko for snitsår på grund af skarpe kanter. Vær forsigtig når du håndterer platforme og andre tilbehørsdele.

FORSIGTIG En sikker fastgørelse af en kolbeholder er kun garanteret i forbindelse med de medfølgende skruer. Sæt de medfølgende skruer i alle kolbeholderens dertil beregnede monteringshuller.

BEMÆRK Hvis skruerne er for lange, kan dette påvirke den udskiftelige platforms stabilitet og funktionalitet. Husk, at du kun må bruge de medfølgende skuer til fastgørelsen.

Kolbeholder

Hver kolbeholder består af en beholder-holdeklemme, en eller to fjedre alt efter kolbeholder-/flaskestørrelse og skruer til fastgørelse på platformen. Anvend udelukkende de skruer, der medfølger til kolbeholderen.



① Fjederbøsning ② Fjeder ③ Skruer ④ Holdeklemme ⑤ Fjeder

Afbildning 81: Kolbeholder med 2 fjedre

- 1. Fastgør evt. fjedrene til holderklemmerne som vist i illustrationerne.
- 2. Fjederbøsningerne skubbes mellem holdeklemmerne som vist i illustrationen. Der anvendes to fjedre til nogle kolbeholdere. Den anden fjeder sættes i omkring den nederste ende af holdeklemmen, efter kolbeholderen er fastgjort til platformen.
- 3. Skru holdeklemmerne fast til platformen med de medfølgende skruer.

<u>Beholder</u>

 Sæt den ønskede beholder forsigtigt i kolbeholderen ved først at trække holdefjederen langt nok væk så beholderens bund kan placeres midt i kolbeholderen. Lad beholderen glide uhindret i den rigtige position og fastgør den til kolbeholderens bredere bund. Fjederen holder beholderens hals sikkert på plads.



 Kontroller, at alle beholdere sidder godt fast, før du tænder apparatet.

Beholderen bør om muligt lukkes med en prop for at forhindre, at stoffer bliver slynget ud under blandingen.

3. 3. 3. Installation af firkantet kolbeholder

FORSIGTIG Biologisk risiko på grund af ødelagte eller utætte beholdere. Forkert installerede tilbehørsdele kan medføre spildte prøver. Kontroller, at tilbehørsdelene er installeret korrekt ved at anvende de rigtige værktøjer og skruer. Til isætningen er en skruetrækker (PH2) med en skaftlængde på 150 mm/6 in nødvendig. (Artikelnummer 75004131). Kontroller, at tilbehørsdelene er placeret korrekt på platformen. Anvend altid en beholder med tilbehør, der er tilpasset den pågældende størrelse. Beholderne skal være intakte og installeret korrekt.

FORSIGTIG Risiko for snitsår på grund af skarpe kanter. Vær forsigtig når du håndterer platforme og andre tilbehørsdele.

Firkantet kolbeholder

Hver firkantet kolbeholder består af en bundplade med monteringshuller og fire fjederbelastede holderklemmer. Der medfølger også skruer til fastgørelse af kolbeholderen på platformen. Anvend udelukkende de skruer, der medfølger til kolbeholderen.



① Monteringshuller ② Fjeder ③ Holderklemme ④ Skruer

Afbildning 82: Firkantet kolbeholder

- 1. Sæt evt. alle fjedre @ i holderklemmerne ③ som vist i Afbildning 82.
- 2. Læg klemmemodulet på platformen og juster monteringshullerne ①.
- 3. Skru klemmemodulet fast på platformen med de medfølgende skruer ④.
- 4. Sæt den firkantede beholder i fra oven.
- 5. Kontroller, at beholderen ligger på bundpladen og sidder godt fast i kolbeholderen.

3. 3. 4. Installation af rørstativ

FORSIGTIG Risiko for snitsår på grund af skarpe kanter. Vær forsigtig når du håndterer platforme og andre tilbehørsdele.

Hvert rørstativ består af to monteringsbøjler (trådramme eller plade) med et fikseringsstifthoved og to fingerskruer pr. rørstativholder til fastgørelse på platformen. Hver rørstativenhed leveres formonteret med et rørstativ og skumindsats(er). Anvend altid de skruer, der medfølger til stativet.



Afbildning 83: Rørstativenhed med monteringsbøjler (trådramme)



- Rørstativ
- Skumindsatser
- ③ Fikseringsstifthoved
- ④ Monteringsholder

Afbildning 84: Rørstativenhed med monteringsbøjler (plade)

- 1. Monter monteringsbøjlerne på platformen ved hjælp af fingerskruerne. Spænd fingerskruerne fast med hånden.
- 2. Sæt rørstativet i monteringsbøjlerne.
- Indstil den ønskede vinkel ved hjælp af fikseringsstifthovederne. Spænd fikseringsstifthovederne fast med hånden.

3. 3. 5. Installation af rørstativholder med justerbar vinkel

FORSIGTIG Risiko for snitsår på grund af skarpe kanter. Vær forsigtig når du håndterer platforme og andre tilbehørsdele.

Rørstativholdere og rørstativer fås i forskellige størrelser. Du finder en liste i afsnittet "1. 2. 5. Rørstativholder med justerbar vinkel" på side 28. Alle rørstativholdere kan indstilles i syv positioner, kan drejes fra 0° i begge retninger og kan låses fast ved 15°, 30° og 45°.

Anvend altid de skruer, der medfølger til stativholderen.



Afbildning 85: Rørstativholder med justerbar vinkel med installeret rørstativ

- 1. Monter stativholderen ① på rysteapparatets platform med de medfølgende skruer
- Spred metalpladerne ③ i begge ender af stativholderens drejeområde ⑥ en smule og sæt rørstativet ② i.
- 3. Sæt rørstativet 2 fast ved forsigtigt at løsne pladerne 3.
- 4. Træk fikseringsstifternes hoveder (5) udad i begge ender af holderen. Stifterne låses op med en 1/4 omdrejning af hovedet.
- Drej drejeområdet

 for stativholderen
 til den ønskede vinkel på 15°, 30° eller 45°.
- 6. Sæt fikseringsstiften (5) i placeringshullet (4) igen og lås stiften fast ved at dreje hovedet 1/4 omdrejning.
- For at fjerne stativet spredes metalpladerne (3) i begge ender af drejeområdet (6), og rørstativet (1) løftes ud.

3. 3. 6. Installation af kolbeholder til mikrotiter-/Deepwell-plader

FORSIGTIG Risiko for snitsår på grund af skarpe kanter. Vær forsigtig når du håndterer platforme og andre tilbehørsdele.



 Deepwell-plade
 Skrue
 Mikrotiterpladens kolbeholder

Afbildning 86: Enheden kolbeholder til mikrotiter-/Deepwell-plader

- 1. Placer mikrotiter-pladerammen på platformen.
- Monter mikrotiter-pladerammen på platformen med de skruer, der medfølger til mikrotiter- / Deepwell-pladesættet.
- 3. Før mikrotiter- eller Deepwell-pladen ind i mikrotiter-pladerammen.
- Kontroller, at mikrotiter- eller Deepwell-pladen sidder fast ved at løfte den forsigtigt.

3. 3. 7. Installation af justerbar beholderholder

FORSIGTIG Risiko for snitsår på grund af skarpe kanter. Vær forsigtig når du håndterer platforme og andre tilbehørsdele.





 ① Monteringshuller ② Justerbare stænger ③ Fikseringsstifthoveder ④ Bakkens sidevæg ⑤ Skridsikker måtte ⑥ Skruer

Afbildning 87: Enhed med enkel højdejusterbar beholderholder

Tilbehør





 ① Monteringshuller ② Justerbare stænger ③ Fikseringsstifthoveder ④ Bakkens sidevæg ⑤ Skridsikker måtte ⑥ Skruer

Afbildning 88: Enhed med dobbelt højdejusterbar beholderholder

Sådan udfører du installationen:

- 1. Løft den skridsikre måtte for at blotlægge monteringshullerne (pos. 1) i Afbildning 88).
- 2. Stil den justerbare beholderholder på platformen og tilpas monteringshullerne.
- 3. Skru den justerbare beholderholder fast på platformen med de medfølgende skruer 6.
- 4. Læg den skridsikre måtte 6 i bunden af den justerbare beholderholder.
- 5. Stil beholderen/beholderne på den skridsikre måtte ⑤.
- 6. Løsn fikseringsstifthovederne 3.
- 7. Indstil de justerbare stængers position [®] således at beholderne er spændt sikkert fast.
- 8. Spænd fikseringsstifthovederne 3 fast igen.
- 9. Kontroller, at beholderen er spændt sikkert fast ved forsigtigt at løfte den.

101 / 141

3. 3. 8. Installation af bægerglasstativer

FORSIGTIG Risiko for snitsår på grund af skarpe kanter. Vær forsigtig når du håndterer platforme og andre tilbehørsdele.



Afbildning 89: Bægerglasstativets enhed

- 1. Læg den nederste plade 2 på platformen.
- 2. Sæt bægerglasstativet ① på den nederste plade.
- 3. Tilpas monteringshullerne efter platformen.
- 4. Skru bægerglasstativets enhed fast med de medfølgende skruer ③.

3. 3. 9. Installation af skilletragtholder

FORSIGTIG Risiko for snitsår på grund af skarpe kanter. Vær forsigtig når du håndterer platforme og andre tilbehørsdele.

BEMÆRK Kontroller, at du anvender de rigtige skruer. Der medfølger to skruetyper. De lange skruer med rød skruesikringslak skal anvendes til sikring af den vertikale stander. De korte skruer uden skruesikringslak må kun anvendes til montering af skilletragtholderen på platformen.

Skilletragtholderen kan monteres på rysteapparatets platform på to grundlæggende måder: horisontalt og vertikalt.



Afbildning 90: Skilletragtholderens monteringspositioner

Den horisontal position ③ i Afbildning 90 kan indstilles i højden for at opnå den skrå position, der er vist i ④. Derudover kan skilletragtholderen kombineres med den vertikale stander (fås som tilbehør) for at få de i ① og ② viste vertikale positioner. Holderen kan også anbringes således, at tragtens hane vender nedad ①, eller omvendt, for at opnå den i ② viste position.

Forberedelse af vertikal montering

Afbildning 91 viser ® skilletragtholder-enheden ® med passende tilbehør til vertikale standere ved siden af hinanden.



Afbildning 91: Skilletragtholder (venstre) med tilbehør til vertikal stander (højre)

- Fastgør skilletragtholderen i den vertikale standers lange ende som vist i ① og ② i Afbildning 90.
- 2. Juster den på følgende måde i forhold til de fire monteringshuller i den vertikale standers tragtholder (@ i Afbildning 91):
 - a. Anvend hullerne til vertikal standard-montering (③ i Afbildning 91) for at montere tragthanen, så den vender nedad.
 - b. Anvend hullerne til vertikal omvendt montering (④ i Afbildning 91) for at montere tragthanen, så den vender opad.
 - c. Skru skilletragtholderen fast på den vertikale stander med de medfølgende skruer (⑤ i Afbildning 91).

Montering på platformen

- 1. Stil skilletragtholder-enheden på platformen.
- 2. Tilpas monteringshullerne efter platformen.
- 3. Skru skilletragtholder-enheden fast på platformen med de skruer, der medfølger til skilletragtholderen (6 i Afbildning 91 og Afbildning 92).

Tilbehør

Installation af tragten



- Skilletragt
- ② Bagerste tragtholder
- ③ Forreste tragtholder

④ Fikseringsstifthoved,

højdeindstilling

⑤ Fikseringsstifthoved,længdeindstilling

6 Skruer uden

skruesikringslak

Afbildning 92: Skilletragtholder-enhed

- Løsn fikseringsstifthovedet til længdeindstilling 6 og skub den forreste tragtholder
 helt fremad.
- 2. Før tragtens bagerste ende ① ind i den bagerste tragtholder ②.
- 4. Indstil den forreste tragtholder ③ i en passende højde og spænd den fast ved at spænde fikseringsstifthovedet til højdeindstilling fast ④. Du kan anbringe skilletragten i en skrå stilling ved at ændre højdeindstillingen.
- 5. Spænd den forreste tragtholder ③ fast ved at spænde fikseringsstifthovedet til længdeindstilling ⑤.

3. 3. 10. Installation af universalbakke

FORSIGTIG Risiko for snitsår på grund af skarpe kanter. Vær forsigtig når du håndterer platforme og andre tilbehørsdele.



Afbildning 93: Universalbakkeenhed

Sådan udfører du installationen:

- 1. Læg universalbakken 2 på platformen.
- 2. Tilpas monteringshullerne efter platformen.
- 3. Skru universalbakken 2 fast med de skruer, der medfølger til universalbakken 3.
- 4. Læg gummimåtten 1 i bakken 2.
- Læg flaskerne på bakken og spænd dem fast med de medfølgende O-ringe ④, se Afbildning 94.

BEMÆRK Det anbefales at bruge O-ringe til omdrejningstal over 100 o/min.



Afbildning 94: Universalbakke med fastspændte beholdere og O-ringe

Tilbehør

BEMÆRK Pas på at O-ringene ikke bliver spændt for meget, når beholderne spændes fast. Når du strækker hver O-ring til monteringen, vil du konstatere et dødt punkt i elasticiteten ud over hvilket, O-ringene ikke kan strækkes mere. Når du har nået dette punkt, bør du ikke forsøge at strække O-ringene mere, ellers går de i stykker.

3. 3. 11. Lægning og anvendelse af klæbemåtten

Klæbemåtter egner sig til beholdere med en bred, flad bund som f.eks. Erlenmeyerkolber, Fernbachkolber, medieflasker, mikrotiterplader, petriskåle, cellekulturkolber, målekolber og bægerglas. Generelt kan alle beholdere, der kan stå oprejst, anvendes, men med et aftagende forhold mellem bundflade og højde reduceres også det maksimalt opnåelige rysteomdrejningstal. Høje beholdere med en lille bund er mindre egnet, se Afbildning 95. Store Erlenmeyerkolber (f.eks. 3000 ml) hæfter mere end mindre (f.eks. 100 ml).



Afbildning 95: Klæbemåtte: Egnede kolbeformer

Der findes generelt to typer klæbemåtter:

- Måtte med dårlig vedhæftning (gennemsigtig) 75004111 og 75004117
 - » nem håndtering ved hjælp af dårligere vedhæftning
 - » særlig velegnet til bægerglas
 - » egnet til alle beholdere op til et maksimalt omdrejningstal på 200 o/min (anvendelse til vævskulturer)
- Måtte med god vedhæftning (sort, gennemsigtig) 75004126 og 75004127
 - » pålidelig beholderfiksering ved hjælp af en bedre vedhæftning
 - » egnet til alle beholdere
 - » maksimalt omdrejningstal på 400 o/min (op til 500 ml-Erlenmeyer)

<u>Værktøj</u>

Den i Afbildning 96 viste applikatorrulle leveres som tilbehør til klæbemåtter. Den er beregnet som et hjælpemiddel til at trykke klæbemåtten fast på platformen med.



Afbildning 96: Applikatorrulle til klæbemåtter

Isætning af tør måtte

- 1. Kontroller, at rysteapparatets platform er ren og tør.
- 2. Fjern beskyttelsesfolien fra den ene side af klæbemåtten.
- 3. Klæb måtten fast på platformen med klæbesiden vendende nedad.
- 4. Fjern beskyttelsesfolien fra den øverste side.

Isætning af fugtig måtte

- 1. Fjern beskyttelsesfolien på begge sider af klæbemåtten.
- 2. Læg måtten i blød i vand eller skyl den grundigt af.
- 3. Lad måtten dryppe af.
- 4. Bred den fugtige måtte ud på den rene platform og placer den som ønsket.
- 5. Lad måtten tørre i 24 timer på platformen før du anvender den.

Bemærkninger til beholderegenskaber

Det maksimale omdrejningstal afhænger at beholdernes type, der skelner mellem størrelse, form, bundform, materiale, overfladeruhed og renhed. Du skal kontrollere disse parametre for at bestemme den optimale måtte til din anvendelse, før du anvender rysteapparatet.

Brug Tabel 53 til at bestemme dit laboratorieudstyrs egnethed til anvendelse med klæbemåtter.
Plast-/glasprodukter	Vurdering af vedhæftningen	Metal	Vurdering af vedhæftningen
Glas (borosilikatglas)	5	Rustfrit stål	5
Polypropylen (PP)	3	Aluminium	5
Polypropylen-Copolymer (PPCO)	3	Elokseret aluminium	4
Polyethylen (HPDE) (LPDE)	3		
Polycarbonat (PC)	4		
Polyethylenterephthalat (PET)	4		
Polyethylenterephthalat- Copolyester (PETG)	4		
Polymethylpenten (PMP)	3		
Polytetrafluorethylen (PTFE)	0*		
Tetrafluorethylen- Hexafluorpropylen- Copolymer (FEP)	0*		
Perfluoralkoxy-Polymerer (PFA)	2		
Silikone	1*		
Vurderinger: 0 = ingen vedha	æftning; 5 = bedste ve	dhæftning/* Anvend i	ikke beholdere med

en vurdering på 0 og 1 (PTFE, FEP og silikone).

 Tabel 53:
 Beholdermaterialemes
 kompatibilitet

BEMÆRK Bemærk venligst, at det maksimale rysteomdrejningstal kun kan opnås med de materialer, der er vurderet til 5 i Tabel 53.

BEMÆRK Ved brug af andre materialer end glas skal du selv teste og bestemme, hvilken type materiale, der er egnet til hvilken/hvilket fyldning, omdrejningstal, funktionstid og temperatur.

<u>Omdrejningstalområde og funktionstidsbegrænsning for glasbeholdere</u> (30 % fyldning)

Grænseværdier for omdrejningstal og funktionstid ved stuetemperatur for glasbeholdere (30 % fyldning). Højere temperaturer reducerer vedhæftningen. Hvis temperaturen ændres, kan beholderne løsne sig fra måtten. Start rysteprocessen med akklimatiserede beholdere.

	Mikrotiter/ DeepWell	Bægerglas			
	96 x 2 ml	25 ml	125 ml	250 ml	500 ml
125 o/min	ingen funktionstidsbeg	grænsning			
150 o/min	72 timer	72 timer	72 timer	72 timer	72 timer
200 o/min	72 timer	72 timer	72 timer	72 timer	72 timer
250 o/min	72 timer				
300 o/min					
350 o/min					
400 o/min					

Tabel 54: Måtte med dårlig vedhæftning - Omdrejningstalområde og funktionstidsbegrænsningved stuetemperatur

	Mikrotiter/ DeepWell	Erlenm	neyer						
	96 x 2 ml	25 ml	125 ml	250 ml	500 ml	1000 ml	2000 ml	3000 ml	5000 ml
125 o/min	ingen funktior	nstidsbeg	grænsnir	ig					
150 o/min	72 timer	72 timer							
200 o/min	72 timer	72 timer							
250 o/min	72 timer	72 timer	72 timer	72 timer	72 timer	72 timer	72 timer	48 timer	24 timer
300 o/min	72 timer	72 timer	72 timer	72 timer	72 timer	72 timer	48 timer	24 timer	
350 o/min	48 timer	48 timer	48 timer	48 timer	48 timer	48 timer	24 timer		
400 o/min	24 timer	24 timer	24 timer	24 timer	24 timer				

Tabel 55: Måtte med god vedhæftning - Omdrejningstalområde og funktionstidsbegrænsningved stuetemperatur

Forberedelse

FORSIGTIG Brug altid beskyttelsesbriller og handsker til håndtering af glasprodukter. Før beholderne placeres på klæbemåtten:

- 1. Kontroller beholderne for mulige skader. Anvend ikke beskadigede beholdere.
- 2. Tør beholderen af med en ren og tør klud for at fjerne alt fugt.
- 3. Fjern fedt og andre urenheder.
- 4. Kontroller, om beholderens bund og klæbemåtten er tør og ren.

FORSIGTIG En næsten usynlig film eller en enkelt vanddråbe på beholderens bund er nok til at medføre, at beholderen mister vedhæftningen og løsner sig fra måtten efter en vis funktionstid.



Afbildning 97: Klæbemåtte: Tør beholderen af

5. Reducer fyldningen eller luk beholderne for at undgå at spilde væsken.

FORSIGTIG Selvom en beholder hæfter perfekt, kan en enkelt dråbe væske, der løber ned af den udvendige side, medføre, at beholderen løsner sig fra måtten efter en bestemt funktionstid.

Isætning af beholdere

- 1. Placer beholderne på måtten og tryk dem nedad.
- 2. Tryk elle træk forsigtigt så hver beholder hæfter godt på måtten.

BEMÆRK Bemærk, at der kan dannes kondensvand ved lave temperaturer eller ved længere tids brug, der muligvis kan medføre, at beholderne løsner sig utilsigtet. Prøve- og beholdertemperaturen bør ikke afvige fra det angivne temperaturområde for rysteapparatet. Du finder yderligere oplysninger om dit rysteapparats mulige temperaturområde i afsnittet "1. 1. Tekniske data" på side 11.

Tilbehør

Fjernelse af beholdere

- 1. Tryk på knappen Stop i displayet eller vent, indtil processen er afsluttet.
- 2. Vent, indtil platformen står helt stille.
- 3. Træk beholderens hals forsigtigt og ensartet til siden og vent et par sekunder. Det kan vare et par sekunder at løsne beholdere med store bundflader.



Afbildning 98: Klæbemåtte: Skub eller træk forsigtigt beholderen af

FORSIGTIG Skub eller træk forsigtigt uden brug af kræfter! Beholderne kan pludseligt løsne sig og spilde indholdet!

4. Hæld en smule vand omkring bunden, hvis en beholder ikke kan løsnes.

Lad måtten lufttørre før du bruger den igen.

Sådan fjernes måtten

1. For at løsne måtten løfter du den i den ene side og trækker den langsomt af platformen som vist i Afbildning 99.



Afbildning 99: Klæbemåtte: Løft og træk måtten væk for at fjerne den

Rengøring og desinfektion

Rengør måtten når vedhæftningen forringes.

FORSIGTIG Anvend ikke andre opløsnings-, vaske- eller rengøringsmidler end et mildt opvaskemiddel uden hudplejende tilsætningsstoffer. Den anbefalede koncentration er maks. 2 ml rengøringsmiddel til 1 liter vand.

Tilbehør

BEMÆRK Hvis måtten kommer i direkte kontakt med koncentreret rengøringsmiddel, reduceres vedhæftningen. I dette tilfælde er det ikke nok at skylle den. Ved direkte anvendelse af rengøringsmidler skal måtten skylles, opblødes i vand i 24 timer og derefter tørre i 24 timer for at gendanne den fulde vedhæftning.

BEMÆRK Desinficer kun med 70 % isopropanol eller 70 % ethanol.

Lettere snavset eller støvet:

- 1. Tør måtterne på platformen fugtigt af fra oven eller desinficer dem.
- 2. Fugt måtten.
- 3. Tør rester af med en svamp.
- 4. Tør måtten fugtigt af med en svamp, der er vædet med rent vand.
- 5. Desinficer evt. måtten.
- 6. Lad måtten tørre i 24 timer.

Meget snavset:

- 1. Fjern måtten fra platformen for at rengøre den fuldstændigt.
- 2. Opblød måtten i vand med fortyndet opvaskemiddel.
- 3. Gnid måtten grundigt af med en svamp for at fjerne alle rester.
- 4. Skyl måtten grundigt af med vand efter rengøringen.
- 5. Desinficer evt. måtten.
- 6. Lad måtten tørre i 24 timer.
- 7. Måtten sættes i som beskrevet i afsnittet "Isætning af tør måtte" på side 108 eller "Isætning af fugtig måtte" på side 108.

3. 3. 12. Installation af gasfordeler

Gasfordeleren giver mulighed for at lede en inert gas som kvælstof eller kuldioxid ind i flere beholdere, når rysteapparatet er i drift (du finder bestillingsoplysninger i afsnittet "1. 2. 12. Generelt tilbehør" på side 42).

Gasfordeleren er en fabriksinstalleret valgmulighed til rysteapparaterne Solaris 2000 I/R og 4000 I/R, der også fås som tilbehør, der skal installeres af kunden, hvis de ikke oprindeligt er installeret på apparatet. Alle rysteapparater af typen Solaris 2000 I/R og 4000 I/R har en kabelgennemføring på en af de to sidevægge (du finder oplysninger i afsnittet "1. 4. Produktoversigt" på side 44) til gennemføring af laboratoriegasforsyningsslangen til fordelerenheden. Gasfordeleren har en indløbsåbning til gastilførslen og otte udløbsåbninger til beholdere.



 Laboratoriegasforsyningsslange

- ② Gasfordelerenhed
- ③ Udløbsåbninger med slanger
- ④ Indløbsåbning
- 5 Fastspændingsskruer

Afbildning 100: Gasfordeler, der er installeret i rysteapparatet Solaris 4000 R

Sådan udfører du installationen:

- 1. Skru gasfordeleren @ fast på bagsiden af rysteapparatets kammer med de medfølgende skruer (5).
- 2. Fjern plastproppen fra den udvendige side af kabelgennemføringstyllen (du finder oplysninger om positionen i afsnittet "1. 4. Produktoversigt" på side 44).
- 3. Skær et kryds i gummityllens membran i rysteapparatets kammer med en hobbykniv.
- 4. Skær et stykke slange til en passende længde til anvendelse som laboratoriegasforsyningsslange ①.

Tilbehør

- 5. Sæt laboratoriegasforsyningsslangen ① fast på krogen på fordelerens indløbsåbning ④.
- 6. Sørg for, at laboratoriegasforsyningsslangen ① ikke når skærmens greb og før den gennem kabelgennemføringens gummitylle og udad.
- 7. Tilslut laboratoriegasforsyningsslangen ① til gasforsyningen.
- Skær 8 slangestykker af og fastgør dem ved fordelerens udløb ③. Anvend Y-forbindelsesstykker ved mere end 8 beholdere og frakobl ubenyttede slanger ved mindre end 8 beholdere for at reducere gastab.
- 9. Anvend egnede propper og tætninger til indføring af slangerne i beholderne.

3.4. Fyldning og tilsigtet anvendelse



Brandfare på grund af udløste kemiske reaktioner.

Start ikke rysteapparatet med omdrejningstal, hvor beholderens indhold kan blive slynget ud.

ADVARSEL

Øg omdrejningstallet langsomt. Før der anvendes kemikalier, anbefaler vi en testkørsel med vand.



Sikkerheden kan blive forringet på grund af forkert fyldning og beskadigede tilbehørsdele.

FORSIGTIG

- Kontroller, at fyldningen (tilbehør og prøver) er fordelt symmetrisk til midten af platformen. Ved brug af en totrins platform skal man være opmærksom på, at den nederste platform skal være fyldt mere end den øverste.
- Overfyld aldrig rysteapparatet. Til maksimal belastning se "Tekniske data" på side 11. Belastningen omfatter vægten af platformen, der er monteret på rysteapparatet, tilbehørsdele og prøver.
- Kontroller, at tilbehørsdelene er installeret korrekt, før du tager rysteapparatet i brug. Følg anvisningerne i afsnittet "Tilbehør" på side 89.



Fare for finger- eller håndskader på grund af nedfaldende låg.

- FORSIGTIG
- Rysteapparatets låg kan lukke utilsigtet og klemme og kvæste fingre og hænder. Dette kan ske, hvis lågets fjeder er defekt, og låget ikke kan holdes i dets helt åbnede position. Brug ikke rysteapparatet med en defekt lågfjeder. Kontakt Thermo Fisher Scientific-kundeservice og få omgående lågfjederen udskiftet.

Påfyldning

Kontroller, at fyldningen inkl. tilbehør og prøver er fordelt symmetrisk omkring midten af platformen. Ved brug af en totrins platform uden udnyttelse af den fuldstændige kapacitet: Der opnås optimale resultater ved fortrinsvis at påfylde den nederste plade og at anbringe fyldningen symmetrisk omkring midten af platformen.

Ved anvendelse af et temperaturstyret rysteapparat skal man sørge for, at fyldningen er placeret således, at der er nok plads til at en kollision og mulig beskadigelse af fyldningen undgås, når låget lukkes, og når det er lukket.



Afbildning 101: Eksempel på en korrekt fyldt platform

_												ā.		i de la compañía de l
•		٠			*		٠	٠		÷	-	•		()
•		*	٠	٠	٠	٠	٠		٠		. "	÷	·(•)	
٠		٠		٠	*	*	*	*	*	٠	*	6	, S	à
			٠			*		*	٠		*		· 20	e(•)}
٠		٠	*		*	*	*	٠	*	•	٠	*	(\cdot)	
٠			٠	*	٠		٠				٠		, N	à
													·	(•)))
						۲				Ø				
	a	*			*		*		•	*	*		K°.	8 W
•		٠		•					٠	٠	٠	۰	. 😴	· A
•		•	*		*		*	*		*	*	*	· ሕ	(•))
•		•	٠		٠	•	٠	٠		٠	٠	٠	-(•)	h 🐨
													100	

Afbildning 102: Eksempel på en forkert fyldt platform

BEMÆRK Den maksimale fyldning omfatter vægten af den platform, der er monteret på rysteapparatet, af tilbehørsdelene og af prøverne. Du finder yderligere oplysninger om den maksimale fyldning i afsnittet "1. 1. Tekniske data" på side 11.

Reglementeret anvendelse

Informationerne om grænseværdierne for omdrejningstal og fyldning er kun vejledende værdier til anvendelse af dette produkt. En rolig drift med få eller ingen vibrationer afhænger af mange faktorer som f.eks. fyldningsvolumen, kolbeholdernes og måtternes type og tilstand, beholdernes placering på platformen, rysteapparatets underlag og det nominelle omdrejningstal.

Hver gang, du klargør en ny eller ukendt kombination eller ændrer en af disse parametre: Øg omdrejningstallet langsomt til 175 o/min og kontroller, om der opstår uønskede svingninger af fyldningen eller om rysteapparatet bevæger sig. Kunden er selv ansvarlig for, at rysteapparatet er sikkert at bruge.

Rysteapparater i Solaris-serien er specielt beregnet til en lav varmeafgivelse og lave vibrationer og egner sig derfor fremragende til brug i en række laboratorieudstyr, herunder klimakamre, inkubatorer og laboratoriekøleskabe. På grund af diverse specifikke omgivelsesbetingelser som bl.a. platformenes og holdernes stabilitet er brugeren selv ansvarlig for, at laboratorieudstyret fungerer sikkert hver gang, det bruges. Øg omdrejningstallet langsomt og vær opmærksom på om apparatet udfører usædvanlige bevægelser.

Sikkert omdrejningstal

Det fyldningsuafhængige sikre omdrejningstal for rysteapparater i Solaris-serien fremgår af Tabel 56. Fyldningen omfatter vægten af den platform, der er monteret på rysteapparatet, af tilbehørsdelene og af prøverne. Du finder yderligere oplysninger om den maksimale fyldning i afsnittet "Tekniske data" på side 11.

Туре	Maks. sikkert omdrejningstal
Solaris 2000	175 o/min
Solaris 4000	175 o/min
Solaris 2000 I	250 o/min
Solaris 2000 R	250 o/min
Solaris 4000 I	250 o/min
Solaris 4000 R	250 o/min

Tabel 56: Maks. sikkert omdrejningstal

Maksimalt omdrejningstal

Ved det højeste omdrejningstal på 525 o/min kan rysteapparater i Solaris-serien bruges med en begrænset fyldning iht. Tabel 57. Fyldningen omfatter vægten af den platform, der er monteret på rysteapparatet, af tilbehørsdelene og af prøverne. Tag højde for de anvendte kolbeholderes og beholderes tilladte omdrejningstalområder, der evt. ikke er kompatible med dette omdrejningstal.

Туре	Maks. sikre fyldninger
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

Tabel 57: Maksimal fyldning ved maksimalt omdrejningstal

BEMÆRK Der skal altid tages højde for, at kolbeholderne begynder at åbne sig med et stigende omdrejningstal og på grund af kraftigere beholderbevægelser sørger for en yderligere forbigående systemsvingning.

Kurver for vægt/omdrejningstal

Følgende kurver for vægt/omdrejningstal hjælper dig med at indstille den optimale kombination af omdrejningstal og fyldning til din tilsigtede anvendelse. Det grønne område viser kombinationerne af omdrejningstal og fyldning, der ikke forårsager nogen eller kun mindre svingninger under rysteprocessen. Det røde område viser kombinationer af omdrejningstal og fyldning, der forårsager kraftige svingninger under rysteprocessen, og som kan medføre en uønsket bevægelse af rysteapparatet. Fordi disse diagrammer kun er en vejledning, skal du passe på, når din kombination af omdrejningstal og fyldning nærmer sig det røde område.

Bemærk, at dine specielle anvendelsesbetingelser kan forårsage, at din fyldning eller dit apparat reagerer på en uønsket måde, før den viste grænselinje er nået. Øg omdrejningstallet trinvist for at undersøge, hvordan din specielle fyldning reagerer. Fyldningen omfatter vægten af den platform, der er monteret på rysteapparatet, af tilbehørsdelene og af prøverne.

BEMÆRK Hver blå horisontal linje i følgende diagrammer angiver en bestemt platforms egenvægt.

Solaris 2000



Afbildning 103: Solaris 2000 - Tilsigtet anvendelse



Solaris 4000

Afbildning 104: Solaris 4000 - Tilsigtet anvendelse

Fyldning og tilsigtet anvendelse



Afbildning 105: Solaris 2000 I / 2000 R - Reglementeret anvendelse



Solaris 4000 I / 4000 R

Afbildning 106: Solaris 4000 I / 4000 R - Reglementeret anvendelse

4. Vedligeholdelse og pleje



ADVARSEL

Fare ved håndtering af farlige stoffer

- Læs og følg Verdenssundhedsorganisationen WHO's Laboratory Biosafety Manual og nationale bestemmelserne, hvis du ryster nogen form for farlige materialer. Hvis der rystes mikrobiologiske prøver af risikogruppe II (iht. Verdenssundhedsorganisationen WHO's Laboratory Biosafety Manual), skal der anvendes aerosoltætte biotætninger. Du finder "Laboratory Biosafety Manual" på Verdenssundhedsorganisationens internetside (www.who.int). For materialer i en højere risikogruppe skal der træffes yderligere sikkerhedsforanstaltninger.
- Hvis giftige eller patogene substanser har kontamineret rysteapparatet eller dele af det, skal du gennemføre passende desinfektionsforanstaltninger ("Dekontaminering" på side 125; "Desinficering" på side 125).
- Afbryd strømforsyningen til rysteapparatet og forlad omgående området, hvis der opstår en farlig situation.



ADVARSEL

Sundhedsskader på grund af smitsomme stoffer

Hvis der utilsigtet kommer væske eller andre materialer under platformen, skal rysteapparatet omgående slukkes, lysnetstikket skal trækkes ud, og platformen skal fjernes ("Platforme" på side 18).

Tør de spildte væsker op iht. de gældende standardprocedurer for laboratorierne. Anvend korrekte personlige værnemidler.

Alle interne indstillinger og reparationer skal udføres af en autoriseret servicetekniker. Rysteapparatets hus må ikke åbnes af brugeren.

Følg alle produktinformationer, der medfølger til det pågældende tilbehør, som indeholder specifikke informationer om korrekt vedligeholdelse og rengøring. Anvend kun følgende oplysninger i dette afsnit som retningslinje.

4.1. Principper

For at beskytte personer, miljøet og materialer er du forpligtet til at rengøre rysteapparatet og tilbehørsdelene regelmæssigt og, om nødvendigt, desinficere det/ dem.

Thermo Fisher Scientific anbefaler, at du rengør og desinficerer dit rysteapparat manuelt mindst én gang om måneden. Den normale rumluft indeholder tusindvis af cirkulerende mikroorganismer, der sætter sig i dit rysteapparat, og som kan udgøre en risiko for dine kulturer.

Thermo Fisher Scientific anbefaler at anvende 70 % ethanol, 70 % isopropanol eller desinfektionsmidler på basis af 10 % eller mindre kvaternært ammonium.

BEMÆRK Mekanismen kan blive beskadiget af indtrængende væsker. Sørg for, at ingen væsker og at især ingen organiske opløsningsmidler kommer i kontakt med mekanismen eller mekanismens leje. Organiske opløsningsmidler nedbryder fedtet i mekanismens leje. Hvis der trænger væsker ind i rysteapparatets dele, der ikke er tilgængelige til rengøring, må rysteapparatet ikke tages i brug. Kontakt kundeservice.

BEMÆRK Ikke tilladte metoder eller midler kan angribe rysteapparatets materialer og medføre funktionsfejl. Brug ikke andre rengørings- eller dekontamineringsmetoder end de metoder, der er beskrevet her, hvis du ikke er sikker på, at de er egnet til materialerne. Anvend kun rengøringsmidler, der ikke beskadiger materialerne. Kontakt producenten af rengøringsmidlet i tvivlstilfælde.

BEMÆRK På temperaturstyrede rysteapparater bør du lade låget stå åbent så rengøringsmiddel og vand kan fordampe.

- Træk lysnetstikket ud før rengøring, desinficering eller dekontaminering.
- Fjern installerede tilbehørsdele og platform(e) fra rysteapparatet før rengøring, desinficering eller dekontaminering.
- Anvend varmt vand med et mildt rengøringsmiddel og en blød klud til rengøring af materialerne. Kontakt Thermo Fisher Scientific i tvivlstilfælde. Skyl derefter af med rent vand og lad apparatet tørre grundigt.
- Anvend aldrig ætsende rengøringsmidler som fosforsyre, klorvand eller skurepulver.
- Anvend kun desinfektionsmiddel med en pH-værdi på 6–8.
- Fjern omgående spildt væske med en fnugfri klud, der er fugtet med et ikkeætsende rengøringsmiddel (se rengøringsmiddelproducentens oplysninger).
- Spildt væske kan løbe ind under platformen. Hvis der kommer spildt indhold under platformen, skal du demontere platformen og rengøre skålen under platformen grundigt. Desinficer eller dekontaminer om nødvendigt.
- Kontroller rysteapparatets komponenter og fjern spildte medier eller snavs.

Kontrol af tilbehøret

BEMÆRK Brug ikke et rysteapparat eller tilbehør med tegn på beskadigelse. Det anbefales at få kontrolleret tilbehørsdelene i forbindelse med en regelmæssig rutinevedligeholdelse for at garantere sikkerheden.

Når tilbehørsdelene er blevet rengjort grundigt, skal de kontrolleres for skader, slitage og korrosion.

Metaldele

Ved tegn på en skade som korrosion, slitage eller revner skal tilbehørsdelene omgående tages ud af drift.

Plastdele

Kontroller disse dele for tegn på ridser, skrammer og revner samt om de er falmet. Hvis en undersøgt del viser tegn på skader, skal den omgående tages ud af drift.

<u>Klæbemåtter</u>

Kontroller klæbemåtterne før hver ibrugtagning for tegn på slitage eller beskadigelse som f.eks. revner, misfarvning, gulning, skørhed, deformering, overfladeslitage eller kemikaliepåvirkninger. Bortskaf omgående klæbemåtter, der viser tegn på slitage eller beskadigelse.

4.2. Rengøring

BEMÆRK Før en rengøringsmetode anvendes, skal brugeren kontakte producenten af rengøringsmidlet for at sikre, at den planlagte metode ikke skader materialerne.

BEMÆRK Mekanismen kan blive beskadiget af indtrængende væsker. Lad ikke væsker og især ikke organiske opløsningsmidler komme i kontakt med mekanismen eller mekanismens leje. Organiske opløsningsmidler nedbryder fedtet i mekanismens leje. Vask apparatets udvendige side med en blød klud og en opløsning af en mild sæbe og vand og skyl det derefter af med rent vand og tør det grundigt.

Du finder yderligere oplysninger om korrekt rengøring af rysteapparatet og de anvendte tilbehørsdele i afsnittet "Principper" på side 123.

BEMÆRK På temperaturstyrede rysteapparater bør du lade låget stå åbent så rengøringsmiddel og vand kan fordampe.

Kontrolpanel

Sådan rengøres betjeningspanelet:

- 1. Træk netstikket ud.
- 2. Rengør kontrolpanelet med en tør mikrofiberklud.
- 3. Fugt om nødvendigt mikrofiberkluden med vand og rengør betjeningspanelet igen.
- 4. Lad alle komponenter lufttørre.

4.3. Desinficering

ADVARSEL Fare ved håndtering af farlige stoffer. Rør ikke ved inficerede dele. Ved berøring af kontaminerede dele er der mulighed for alvorlige infektioner. Infektiøst materiale kan trænge ind i rysteapparatet, hvis beholderen går i stykker, eller hvis væsken spildes. Kontroller, i tilfælde af en kontaminering, at ingen er udsat for fare. Desinficer omgående de pågældende dele.

BEMÆRK Materialer kan blive beskadiget på grund af uegnede desinficeringsmetoder eller -midler. Kontroller, at desinfektionsmetoden eller -midlet ikke skader materialerne. Kontakt producenten af desinfektionsmidlet i tvivlstilfælde. Følg sikkerhedsanvisningerne og brugsanvisningerne til de anvendte desinfektionsmidlet.

- 1. Tør alle komponenter og flader af med 70 % ethanol alt efter den nødvendige desinfektionsintensitet. Fugt ikke områder med blotlagte elektronikkomponenter.
- 2. Lad alle komponenter lufttørre.

BEMÆRK På temperaturstyrede rysteapparater bør du lade låget stå åbent så rengøringsmiddel og vand kan fordampe.

Du er selv ansvarlig for, at den dekontamineringsgrad, der opfylder dine krav, opnås.

4.4. Dekontaminering

ADVARSEL Fare ved håndtering af farlige stoffer. Rør ikke ved kontaminerede dele. Ved berøring af kontaminerede dele er der mulighed for kontaminering. Kontamineret materiale kan trænge ind i rysteapparatet, hvis beholderen går i stykker, eller hvis væsken spildes. Kontroller i tilfælde af en kontaminering, at ingen er udsat for fare. Dekontaminer omgående de berørte dele.

BEMÆRK Materialer kan blive beskadiget på grund af uegnede dekontamineringsmetodereller midler. Kontroller, at dekontamineringsmetoden eller -midlet ikke skader materialerne. Kontakt producenten af rengøringsmidlet i tvivlstilfælde. Følg sikkerhedsanvisningerne og brugsanvisningerne til de anvendte dekontamineringsmidlet.

Thermo Fisher Scientific anbefaler følgende metode.

- 1. Tør alle komponenter og flader af med 70 % ethanol. Fugt ikke områder med eksponeret elektronik.
- 2. Lad alle komponenter lufttørre.

BEMÆRK På temperaturstyrede rysteapparater bør du lade låget stå åbent så rengøringsmiddel og vand kan fordampe.

Du er selv ansvarlig for, at den dekontamineringsgrad, der opfylder dine krav, opnås.

4.5. Autoklavering

BEMÆRK Overskrid aldrig de tilladte værdier for autoklaveringstemperatur og -varighed. Kemiske tilsætningsstoffer i dampen er ikke tilladt.

Før autoklaveringen:

- Demonter alle tilbehørsdele.
- Skyl alle spor af kemikalier eller rengøringsmidler grundigt af. Gennemfør en afsluttende skylning med destilleret vand.

Du finder informationer om autoklaverbarhed i følgende tabel:

Artikel	Autoklaverbar	Tekniske data
Rysteapparat	Nej	-
Platform	Ja	121 °C, 20 min.
Kolbeholder (uden fjederbøsninger)	Ja	121 °C, 20 min.
Skruetrækker	Nej	-
Kolbeholder til mikrotiter-/Deepwell-plader	Ja	121 °C, 20 min.
Klæbemåtte	Nej	-
Bægerglasstativ	Ja	121 °C, 20 min.
Skilletragtholder	Ja	121 °C, 20 min.
Rørstativ – Metaldele	Ja	121 °C, 20 min.
Rørstativ – Plastdele	Nej	-
Rørstativ – Skumindsatser	Nej	-
Justerbar beholderholder – Metaldele	Ja	121 °C, 20 min.
Justerbar beholderholder – Skridsikker måtte	Nej	-
Universalbakke – Metalbakke	Ja	121 °C, 20 min.
Universalbakke – Gummimåtte	Nej	-
Rørstativholder med justerbar vinkel – Metaldele	Ja	121 °C, 20 min.
Rørstativholder med justerbar vinkel – Plastdele	Nej	-
Rørstativholder med justerbar vinkel – Skumindsatser	Nej	-

 Tabel 58: Materialers autoklaverbarhed

Kontroller, at den nødvendige sterilitet i overensstemmelse med dine egne krav er opnået.

Udseendet og farven kan ændre sig en smule efter autoklaveringen.

4.6. Temperaturkalibrering

Et temperaturstyret rysteapparats nominelle temperaturværdi kan kalibreres med en temperaturmåler med den nødvendige målenøjagtighed for at sikre en fuldstændig reproducerbarhed af den inkuberede (kun opvarmede) eller kølede (kølede og opvarmede) rysteproces.

4. 6. 1. Pålidelige metoder til temperaturkalibrering

Følg anvisningerne nøje før og under temperaturkalibreringen af rysteapparatet. Enhver forstyrrelse eller afvigelse fra den anbefalede pålidelige metode kan medføre, at prøvetemperaturen ikke stabiliseres, og at kalibreringen mislykkes.

Pålidelige metoder:

- Anvend en kolbe med en kapacitet på 250 ml.
- Fyld kolben med nøjagtigt 100 ml vand.
- Luk låget fuldstændigt og hold det lukket under hele kalibreringen.
- Vent tilstrækkelig længe for at sikre at prøvetemperaturen er fuldstændig stabiliseret.
- Hold øje med at omgivelsesbetingelserne som temperatur og luftfugtighed ikke ændrer sig under kalibreringen. Selv små ændringer som at åbne en dør eller et vindue eller at ændre klimaanlæggets indstillinger kan påvirke resultatet af den igangværende kalibreringsproces.
- Det anbefales at kalibrere rysteapparatet en gang årligt.

4. 6. 2. Gennemførelse af temperaturkalibrering

Sådan gennemfører du kalibreringen:

- 1. Kalibreringen kan startes på tre måder:
 - » Ved afslutningen af den i "Grundindstilling" på side 54 beskrevne grundindstilling har du mulighed for at gennemføre temperaturkalibreringen i stedet for at afslutte grundindstillingen.
 - » Rysteapparatet viser en påmindelse om den årlige kalibrering.
 - » Du kan derefter starte kalibreringen ved at trykke på symbolet Indstillinger på navigationslinjen og derefter på knappen Kalibrering længere nede på det viste skærmbillede Temp.-kalibrering.
 - » Du kan trykke på symbolet **Indstillinger** på navigationslinjen og derefter på knappen **Kalibrering** for at starte processen, når det er nødvendigt.

BEMÆRK Kalibreringspåmindelsen kan deaktiveres ved at klikke på afkrydsningsfeltet Deaktiver kalibreringsmeldinger på siden Indstillinger. 2. Vent, indtil det i Temp.-kalibrering viste skærmbillede Afbildning 107 vises.



Afbildning 107: Kalibrering: Startskærmbillede

3. Tryk på knappen Start kalibrering for at starte kalibreringen.



Afbildning 108: Kalibrering: Opfordring til fyldning af en Erlenmeyerkolbe

- 4. Følg anvisningerne på skærmen: Fyld en 250 ml Erlenmeyerkolbe med 100 ml vand.
- 5. Sæt kolben i en kolbeholder og fastgør kolbeholderen ca. midt på rysteapparatets platform.
- Hæng en temperaturmålers føler med den nødvendige målenøjagtighed i kolben.
 Føleren bør være nedsænket i væsken, men ikke komme i kontakt med kolbens bund eller sider.
- 7. Luk kolben for at forhindre at væsken spildes.
- 8. Luk låget over rysteapparatets platform.

9. Tryk på knappen Næste for at fortsætte.



Afbildning 109: Kalibrering: Opfordring til at starte kalibreringen

BEMÆRK Rysteapparatet starter omgående rysteprocessen, så snart du trykker på knappen Start.

10. Følg anvisningerne på skærmen: Tryk på knappen Start for at anvende den forudindstillede temperatur eller tryk på feltet Nominel temperaturværdi og indtast en anden ønsket temperatur før du trykker på knappen Start.

Skærmbilledet viser, hvordan rysteapparatets kammer bliver opvarmet (eller afkølet) til den valgte kalibreringstemperatur.



Afbildning 110: Kalibrering: Stabilisering af kammeret til kalibreringstemperatur

11. Vent mindst to og en halv time (150 minutter), indtil kammertemperaturmålingen har nået en stabil tilstand.

En stabil tilstand vises med meldingen **Temperaturen i prøverummet er stabiliseret** som vist i Afbildning 111 nedenfor.



Afbildning 111: Kalibrering: Kammer stabiliseret til kalibreringstemperatur

- 12. Følg anvisningerne på skærmen:
 - » Sammenlign regelmæssigt temperaturvisningen på dit termoelement så snart det skærmbillede, der er vist i Afbildning 111, vises. Når du mener, at temperaturen er stabil, kan du enten trykke på knappen **Rediger** for at korrigere den eller på knappen **Afslut** for at afslutte kalibreringsprocessen.



Afbildning 112: Kalibrering: Korrektion af rysteapparatets temperaturvisning

- 13. Korriger den af rysteapparatet viste temperatur til måleudstyrets værdi ved hjælp af det viste indstillingshjul.
- 14. Tryk på knappen Gem for at gemme dine ændringer.

15. Meldingen **Temperaturkalibrering er afsluttet** viser, at kalibreringen blev afsluttet uden fejl.



Afbildning 113: Kalibrering: Temperaturkalibrering afsluttet uden fejl

Du kan trykke på knappen Kalibrer igen for at gentage kalibreringsprocessen.

4. 6. 3. Temperaturkalibreringsforespørgsel

Du kan til enhver tid hente kalibreringstemperaturen:

- Tryk på navigationslinjen på symbolet Indstillinger og derefter på knappen Kalibrering længere nede på det viste skærmbillede Temp.-kalibrering.
- 2. Det i Afbildning 107 viste skærmbillede **Temp.-kalibrering** vises. Dette skærmbillede viser den aktuelle **Kalibrerede rysteapparat-temperatur**.

BEMÆRK På temperaturstyrede rysteapparater kan du alternativt aflæse kalibreringstemperaturen på skærmen, hvor du indstiller temperaturen. Et eksempel er vist i afsnittet "Indstilling af temperaturen" på side 65.

4.7. Firmwareinstallation

Hvis du trykker på knappen Firmware i skærmbilledet Filer og info (se "Filer og info" på side 78), bliver du opfordret til at indtaste administrator-adgangskoden, hvorefter det i Afbildning 114 viste skærmbillede Firmware vises. På dette skærmbillede kan du installere ny firmware fra en USB-nøgle.

Alt efter serviceaftalen får du muligvis et download-link til ny firmware fra din servicetekniker. Sørg for kun at installere firmware fra autoriserede kilder. Ved brug af software fra ikke-autoriserede kilder kan Thermofisher Scientific's garanti bortfalde.

- 1. Download ZIP-filen ved at klikke på linket som du har fået fra din servicetekniker.
- 2. Pak ZIP-filen ud på en tom USB-nøgle.

BEMÆRK Kontroller, at der er nok lagerplads på USB-nøglen til at ZIP-filen kan pakkes ud.

 Hvis rysteapparatet stadig kører, stopper du det ved at trykke på knappen Stop i displayet. 4. Kvitter for alle evt. alarmmeldinger som beskrevet i afsnittet "Driftstilstand" på side 67.

BEMÆRK Hvis du forsøger at gennemføre en opgradering, mens rysteapparatet kører, får du følgende melding: **En opgradering kan ikke gennemføres, mens shakeren er i drift.** Denne melding indeholder en knap **Stop shaker**, med hvilken du kan afslutte processen og gennemføre opgraderingen.

- 5. Tryk på navigationslinjen på symbolet **Filer og info**.
- 6. Rul helt ned på skærmbilledet Filer og info for at få vist knappen Firmware.
- 7. Tryk på knappen **Firmware** for at åbne skærmbilledet Firmware, der viser de aktuelt installerede firmwareversioner.
- 8. Sæt USB-nøglen i en af rysteapparatets USB-indgange. USB-indgangenes placering fremgår af afsnittet "Produktoversigt" på side 44.
- 9. Indtast administrator-adgangskoden.



Afbildning 114: Skærmbilledet Firmware

10. Tryk på knappen Gennemfør firmwareopgradering.

BEMÆRK Hvis rysteapparatet ikke finder en USB-nøgle, bliver du opfordret til følgende: **Sæt en USB-nøgle med den firmware, der skal installeres, i for at starte opgraderingen**. Sæt nu USB-nøglen i eller kontroller, om den er sat rigtigt i USB-indgangen og afhjælp problemet.

11. Hvis en firmwareopgradering skal gennemføres, vises den i Afbildning 115 viste liste på skærmbilledet Firmwareopgradering. Tryk på firmware-elementet for hvilket du vil gennemføre en opgradering.

	^	Firmware Upgrade
Select firmware to upgrade:		Select firmware to upgrade:
O H.M.I.		C H.M.I.
Main Controller		Main Controller
O Parameters		O Parameters
Select		Select

Afbildning 115: Valg af firmware-element, der skal installeres

BEMÆRK Firmware-elementer i den sidste nye version er grå og kan ikke vælges. Hvis der ikke er nogen elementer at opdatere, får du meldingen **Der er i øjeblikket ingen nye opgraderinger til softwaren**.

- 12. Tryk på knappen Vælg.
- Hvis USB-nøglen indeholder mere end én firmwareversion, bliver du opfordret til at vælge en. Tryk på den ønskede firmwareversion og derefter på knappen Vælg.



Afbildning 116: Valg af firmwareversion til installationen

BEMÆRK Hvis du er usikker på, hvilken version du skal installere, bedes du kontakte serviceteknikeren, der har stillet download-linket til rådighed.

14. Der vises en liste over de elementer, som der skal gennemføres en opgradering af. Tryk på knappen **OK** for at starte opgraderingen.



Afbildning 117: Liste over de elementer, som der skal gennemføres en opgradering af

BEMÆRK Brugergrænsefladen administrerer den gensidige afhængighed for dig: Du vælger et firmware-element til opdateringen, og softwaren tilføjer automatisk alle nødvendige elementer til installationen.

BEMÆRK Hvis du vælger valgmuligheden H.M.I. fra den i Afbildning 115 viste liste for kun at opdatere GUI-firmwaren, reagerer GUI-displayet i op til et minut efter starten af firmwareopgraderingen ikke længere på, at der trykkes på brugergrænsefladens aktive område. Sluk <u>ikke</u> for rysteapparatet i dette tidsrum, ellers kan opgraderingen ikke afsluttes uden fejl. Rysteapparatet starter igen efter kort tid og er derefter klar til brug igen.

BEMÆRK Hvis du vælger to eller tre valgmuligheder fra den liste, der vises i Afbildning 115 for at gennemføre en opgradering af <u>flere</u> firmware-komponenter, må rysteapparatet på intet tidspunkt være slukket under hele opgraderingen.

15. Der vises forskellige meldinger, der informerer dig om installationens status som vist i Afbildning 118.



Afbildning 118: Meldinger om installationen af firmwareopgraderingen

16. Hvis du bliver opfordret til at genstarte rysteapparatet, skal du slukke for og tænde apparatet igen med tænd-/sluk-kontakten.

4.8. Udskiftning af platformen

Du finder den grundlæggende fremgangsmåde til montering af en platform i afsnittet "3. 3. 1. Montering af platformen" på side 90. Du finder en fuldstændig liste over reserve-platforme til hvert rysteapparat i afsnittet "1. 2. 1. Platforme" på side 18.

BEMÆRK Temperaturstyrede rysteapparater må kun anvendes med de platforme, der er forudinstalleret på fabrikken.

4.9. Service

Thermo Fisher Scientific anbefaler, at rysteapparatet og tilbehøret vedligeholdes af en autoriseret servicetekniker én gang årligt. Serviceteknikeren kontrollerer følgende:

- Elektriske installationer
- Opstillingsstedets egnethed
- Sikkerhedssystem
- Anvendte tilbehørsdele
- Kolbeholdernes, platformenes og andre tilbehørsdeles fastgørelse på rysteapparatet

Før service bør rysteapparatet og tilbehørsdelene være rengjort og dekontamineret grundigt for at sikre en fuldstændig og sikker inspektion.

Thermo Fisher Scientific tilbyder inspektions- og serviceaftaler i forbindelse med disse eftersyn. Eventuelle nødvendige reparationer udføres gratis inden for garantiperioden og mod betaling, når garantien er udløbet. Dette er kun gældende, hvis det udelukkende er serviceteknikere fra Thermo Fisher Scientific, der har foretaget indgreb i rysteapparatet.

4.10. Forsendelse og bortskaffelse

ADVARSEL Sundhedsskader på grund af smitsomme stoffer. Når du tager rysteapparatet og tilbehørsdelene ud af drift for at bortskaffe det/dem, skal du rengøre dem og evt. desinficere eller dekontaminere dem. Kontakt Thermo Fisher Scientific kundeservice i tvivlstilfælde.

Overhold nationale bestemmelser om bortskaffelse af rysteapparatet. Kontakt Thermo Fisher Scientific-kundeservice for at bortskaffe rysteapparatet. Du finder kontaktoplysningerne på bagsiden af denne vejledning, eller på <u>www.thermofisher.</u> <u>com</u>.

Læs og følg oplysningerne om transport og forsendelse ("Transport" på side 52 og "Forsendelse" på side 60).

5. Årsager til fejl og fejlafhjælpning

Nr.	Betegnelse	Løsninger	Symbol
Fejl			
11999	Der er opstået en intern fejl.	Der er opstået en intern fejl. Genstart rysteapparatet ved at slukke for det og tænde det igen med tænd-/sluk- kontakten. Kontakt en servicetekniker, hvis fejlmeldingen fortsat vises.	
Advarsle	r		
3	Temperaturkalibrering skal udføres. Der er gået 1 år.	Den etårige gyldighedsperiode er udløbet. Forny temperaturkalibreringen som beskrevet i afsnittet "4. 6. Temperaturkalibrering" på side 127.	CAL
4	Strømforsyning genetableret og automatisk genstart aktiveret.	Der er opstået et strømsvigt under det sidste forløb. Efter genetablering af strømforsyningen blev forløbet automatisk genoptaget.	Ŧ
6	Sensorerne for omgivelsestemperatur har svigtet.	Temperaturen, der er målt af omgivelsestemperatursensoren, er ikke plausibel.	III
10	Omdrejningstal for ventilator 1 over grænseværdien.	Omdrejningstallet for ventilator 1 er over grænseværdien.	•2
11	Omdrejningstal for ventilator 2 over grænseværdien.	Omdrejningstallet for ventilator 2 er over grænseværdien.	•2
12	Omdrejningstal for ventilator 3 over grænseværdien.	Omdrejningstallet for ventilator 3 er over grænseværdien.	•2
13	Omdrejningstal for ventilator 4 over grænseværdien.	Omdrejningstallet for ventilator 4 er over grænseværdien.	•2
20	Låget har været åbnet længe.	Luk rysteapparatets låg for at bibeholde kammertemperaturen.	4
Alarmme	ldinger		
1	Alarm overtemperatur	Kammertemperaturen har overskredet den øvre alarmværdi. Kontroller dine prøver, omgivelsesbetingelser og/eller rediger dine indstillinger.	
2	Alarm undertemperatur	Kammertemperaturen har overskredet den nedre alarmværdi. Kontroller dine prøver, omgivelsesbetingelser og/eller rediger dine indstillinger.	
3	Den automatiske genstart efter et strømsvigt mislykkedes.	Der er opstået et strømsvigt under det sidste forløb. Den automatiske genstart kunne ikke gennemføres.	Ċ
5	Drevacceleration for langsom. Det ønskede nominelle omdrejningstal kan ikke nås.	Det ønskede nominelle omdrejningstal kunne ikke nås rettidigt. Kontroller dine indstillinger og/eller platformens fyldning (tilbehør og prøver).	\odot

Forsendelse og bortskaffelse

Nr.	Betegnelse	Løsninger	Symbol
7	Målingen af omdrejningstallet på ventilator 1 melder en uventet tilstand.	Sluk apparatet og tænd det igen. Kontakt en servicetekniker, hvis fejlmeldingen fortsat vises.	2
8	Målingen af omdrejningstallet på ventilator 2 melder en uventet tilstand.	Sluk apparatet og tænd det igen. Kontakt en servicetekniker, hvis fejlmeldingen fortsat vises.	2
9	Målingen af omdrejningstallet på ventilator 3 melder en uventet tilstand.	Sluk apparatet og tænd det igen. Kontakt en servicetekniker, hvis fejlmeldingen fortsat vises.	•
10	Målingen af omdrejningstallet på ventilator 4 melder en uventet tilstand.	Sluk apparatet og tænd det igen. Kontakt en servicetekniker, hvis fejlmeldingen fortsat vises.	•
12	Fejl ved drevstart - intet hastighedssignal modtaget.	Rysteapparatets platform er blokeret. Kontroller, at der er tilstrækkelig plads omkring platformen og/eller reducer fyldningen (tilbehør og prøver) på platformen. Tryk derefter på START igen på rysteapparatet. Kontakt en servicetekniker, hvis fejlmeldingen fortsat vises.	
24	Fejl ved måling af omdrejningstal under et forløb.	Usædvanlig ændring af omdrejningstal konstateret. Kontroller fyldningen (tilbehør og prøver) og/eller kolbeholderne på rysteapparatets platform. Tryk derefter på START igen på rysteapparatet. Kontakt en servicetekniker, hvis fejlmeldingen fortsat vises.	
26	Målingen af omdrejningstallet har konstateret en uventet stilstand under et forløb.	Kontroller, at der er tilstrækkelig plads omkring platformen, og/eller reducer fyldningen (tilbehør og prøver) på platformen. Kontroller, at sikringen på bagsiden af rysteapparatet ikke har slået fra ("Sikring" på side 47). Tryk derefter på START igen på rysteapparatet. Kontakt en servicetekniker, hvis fejlmeldingen fortsat vises.	٢
82	Motorstrømsmålingen har konstateret en overbelastning.	Motor-overstrøm konstateret. Påfyld eller tøm ikke platformen mens apparatet kører. Reducer omdrejningstallet eller tilpas fyldningen (tilbehør og prøver) på platformen.	— 0
83	Motorstrømmens måleværdier er uden for det tilladte område.	Motor-overstrøm konstateret. Påfyld eller tøm ikke platformen mens apparatet kører. Reducer omdrejningstallet eller tilpas fyldningen (tilbehør og prøver) på platformen.	0

Tabel 59: Liste over fejlmeldinger, advarsler og alarmmeldinger

BEMÆRK Hvis der vises en fejlmelding, der ikke findes i denne tabel, bedes du henvende dig til en servicetekniker.

GPL (General Public License, offentlig licens)

En del af apparatets software anvender open source-software, der er udgivet under GPL, LGPL eller en anden open source-licens. Mere specifikt drejer det sig om de libraries, der er anført i tabellen. Kildekoden til de anvendte libraries (libraries fra tredjepartsudbydere) kan købes hos Thermo Fisher Scientific, hvis dette er angivet i den pågældende licens. De pågældende licensbetingelser for den anvendte open source-software er den del af den kildekode-pakke, der er stillet til rådighed.

Library	Version	Anvendelse via	Licens
Qt	5.8	BSP	LGPLv3
Log4Cplus	1.2.0	Арр	Apache Public- licens v2 / BSD- licens med to klausuler
boost	1.72.0	Арр	Boost-licens 1.0
json11	1.0.0	Арр	MIT-licens

Libraries fra tredjepartsudbydere

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

Indeks

Α

Advarsler 70 Alarmmeldinger 70 Apparatbetegnelse 77 Autoklavering 126 Automatisk genstart 74 Årsager til fejl og fejlafhjælpning 136

В

Betjening 61 Bortskaffelse 135

D

Dekontaminering 125 Den lydløse tilstands varighed 71 Desinficering 125 Diagrammer 86 Direktiver 43 Display 74 Driftsparametre 74 Dvaletilstand 76

Е

Ethernet 47

F

Filer og info 78 Forsendelse 60, 135 Fyldning 116

G

GPL (General Public License, offentlig licens) 138 Grafisk brugergrænseflade 61 Grundindstilling 54

Η

Hændelseslog 85

I

Indstillinger 69 Indstillingsværdi for omdrejningstal 74 Installation af bægerglasstativer 102 Installation af firkantet kolbeholder 95 Installation af gasfordeler 114 Installation af justerbar beholderholder 100 Installation af kolbeholder til mikrotiter-/Deepwell-plader 99 Installation af rørstativ 97 Installation af rørstativ 97 Installation af rørstativholder med justerbar vinkel 98 Installation af skilletragtholder 103 Installation af universalbakke 106 Isætning af kolbeholder og beholder 93

Κ

Kontrol af tilbehøret 124

L

Leveringsomfang 49 Lydstyrke 70 Lysnettilslutning 47 Lysstyrke 74 Lægning og anvendelse af klæbemåtten 107

Μ

Maksimalt omdrejningstal 119 Metaldele 124 Montering af platformen 90

Ν

Nominel temperaturværdi 74

0

Opstillingssted 50

Vedligeholdelse 122

Ρ

Plastdele 124 Platforme 18 Pleje 122 Produktoversigt 44 Programmer 79 Påfyldning 117

R

Region 77 Reglementeret anvendelse 5, 118 Rengøring 124

S

Signalord og symboler 6 Sikkert omdrejningstal 118 Sikring 47 Sprog 75 Standarder 43 Symboler, der anvendes i brugsanvisning 7

Т

Tekniske data 11 Tekniske specifikationer 11 Temperaturkalibrering 127 Tid 76 Tidstilstand 74 Tilbehør 17, 89 Tilsigtet anvendelse 116 Tilslutninger 46 Totrins platforme 92 Transport 52

U

Universalplatforme 91 USB 48

V

thermo scientific





Fremstillet til

Thermo Electron LED GmbH Zweigniederlassung Osterode Am Kalkberg, 37520 Osterode am Harz Germany Oprindelsesland: 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 er den originale brugsanvisning. Denne brugsanvisning er en oversættelse af den originale brugsanvisning.

thermofisher.com/shaker

© 2019–2023 Thermo Fisher Scientific Inc. Alle rettigheder forbeholdes.

Hvis intet andet udtrykkeligt er beskrevet, tilhører alle varemærker Thermo Fisher Scientific Inc. og deres datterselskaber. Ikke alle produkter fås i alle lande. Kontakt din lokale forhandler for yderligere oplysninger.

De billeder, der er offentliggjort i denne vejledning, er kun beregnet som reference. Indstillinger og sprog, der er vist i billederne, kan afvige fra Deres centrifuge.

Australien +61 39757 4300

Østrig +43 1 801 40 0

Belgien +32 53 73 42 41

Kina +800 810 5118 **eller** +400 650 5118

Frankrig +33 2 2803 2180

Tyskland nationalt, gratis 0800 1 536 376

Tyskland internationalt +49 6184 90 6000 Indien +91 22 6716 2200

Italien +39 02 95059 552

Japan +81 3 5826 1616

Holland +31 76 579 55 55

New Zealand +64 9 980 6700

Nordiske lande/Baltikum/ SNG-lande +358 10 329 2200

Rusland +7 812 703 42 15 **Spanien/Portugal** +34 93 223 09 18

Schweiz +41 44 454 12 12

Storbritannien/Irland +44 870 609 9203

USA/Canada +1 866 984 3766

Andre asiatiske stater +852 2885 4613

Andre lande +49 6184 90 6000

