

# Thermo Scientific Solaris skakapparater

## Bruksanvisningar

70900197-е • 2024-11



# Innehåll

Förord		5
Om bruks	sanvisningen	5
Korrekt a	nvändning	5
	Öppen skakapparat	5
	Inkuberad skakapparat (uppvärmd) och kyld skakapparat	
	(kyld och uppvärmd)	5
	Kunskapskrav	5
Signalord	och symboler	6
	Symboler som används på skakapparaten och dess	~
		ю 7
		-
Sakerhet	sinstruktioner	7
1. Specifik	ationer	1
1.1. Tek	niska data	1
	Thermo Scientific Solaris 2000 skakapparat1	1
	Thermo Scientific Solaris 4000 Skakapparat1	2
	Thermo Scientific Solaris 2000 I skakapparat1	3
	Thermo Scientific Solaris 4000 I skakapparat1	4
	Thermo Scientific Solaris 2000 R skakapparat1	5
	Thermo Scientific Solaris 4000 R skakapparat1	6
1. 2. Till	pehör 1	7
1. 2. 1.	Plattformar 1	8
1. 2. 2.	Kolvhållare	9
1. 2. 3.	Hållare för mikrotiterplatta/deepwellplatta 2	23
1. 2. 4.	Provrörsställ	24
1. 2. 5.	Provrörsställhållare med justerbar vinkel	28
1. 2. 6.	Justerbara kärlhållare	30
1.2.7.	Kvadratiska mediahållare	32
1. 2. 8.		34
1.2.9.		58
1.2.10	Klistormattor	າສ 10
1 2 10		12
1. 2. 12		r 🗠

1. 3. Dire	ektiv och standarder	43
1.4. Pro	duktöversikt	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.	Anslutningar	46
2. Transpo	ort och uppställning	49
2.1. Upp	backning	49
	Leveransomfattning	49
2. 2. Upp	oställningsplats	50
2. 3. Trai	nsportera	52
	- Hantering	52
2.4. Avv	ägning	53
2. 5. Nät	anslutning	53
2.6. För	sta start	54
2.7. För	varing	59
2.8 Lov	orans	50
2.0. Lev		55
3. Drift		60
3. 1. Strö	òm på/av	60
3. 2. Gra	fiskt användargränssnitt	60
3. 2. 1.	Grundläggande driftinställningar	61
3. 2. 2.	Status	66
3. 2. 3.	Inställningar	68
3. 2. 4.	Program	78
3. 3. Till	oehör	88
3. 3. 1.	Montera plattform	89
3. 3. 2.	Montera kolvhållare och kärl	92
3. 3. 3.	Montera kvadratiska hållare	94
3. 3. 4.	Montera provrörsställ	96
3. 3. 5.	Montera provrörsställhållare med justerbar vinkel	97
3. 3. 6.	Montera hållare för mikrotiterplatta/deepwellplatta	98
3. 3. 7.	Montera justerbar kärlhållare	99

3 3 8 Montera bägarställ	101
3 3 9 Montera separertratthållare	102
3 3 10 Montera universalbricka	102
2 2 11 Montera och använda klistormatta	100
3. 3. 12. Montera gosfördelare	112
	115
3. 4. Lastning och normal anvandning	115
Lastning	116
Normal användning	117
4. Underhåll och skötsel	121
4. 1. Allmän information	122
Inspektera tillbehör	123
4. 2. Rengöring	123
Pekskärm	123
4. 3. Desinficering	124
4.4. Dekontaminering	124
4. 5. Autoklavering	125
4. 6. Temperaturkalibrering	126
4. 6. 1. Etablerad praxis för temperaturkalibrering	126
4. 6. 2. Procedur för temperaturkalibrering	126
4. 6. 3. Temperaturkalibreringsvärden	130
4. 7. Installera inbyggd programvara	130
4. 8. Plattformsbyte	134
4. 9. Service	134
4. 10. Leverans och avyttring	134
5. Felsökning	135
GPL (General Public License)	137
	120
	130

## Förord

## Om bruksanvisningen

Den här bruksanvisningen innehåller beskrivningar av följande skakapparater och kompatibla tillbehör ("1. 2. Tillbehör" på sidan 17):

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

Bruksanvisningen innehåller inte beskrivningar av de olika produkterna utifrån deras fullständiga produktnamn, utan baseras istället på olika klasser av skakapparater:

- Öppen skakapparat
- Inkuberad skakapparat (uppvärmd)
- Kyld skakapparat (kyld och uppvärmd)

## Korrekt användning

## Öppen skakapparat

En enhet som används inom forskning och tillverkning för att skaka lösningar i en 2D-omkretsbana och samtidigt styra hastighet och tid i tillämpningar.

## Inkuberad skakapparat (uppvärmd) och kyld skakapparat (kyld

#### och uppvärmd)

Utrustning som används inom forskning och tillverkning för att skaka lösningar i en 2D-omloppsbana och samtidigt styra hastighet, tid och temperatur i tillämpningar.

#### Kunskapskrav

Skakapparaten får endast användas av utbildad personal.

## Signalord och symboler

Signalord	Risknivå
VARNING	Avser en potentiellt farlig situation, vilken kan leda till dödsolycka eller allvarlig personskada om den inte undviks.
SE UPP	Avser en potentiellt farlig situation, vilken kan leda till smärre eller moderat personskada om den inte undviks.
OBS	Avser information som är viktig, men inte riskrelaterad.

## Symboler som används på skakapparaten och dess tillbehör

Beakta informationen i denna bruksanvisning för att hålla dig själv och din miljö säker.

<b>(</b>	Se bruksanvisningen
	Koppla ur stickkontakten
	Allmän fara
	Fara orsakad av varm yta
	Risk för finger- eller handskador
	orsakade av mekaniska delar som
	oavsiktligt stängs.

## Symboler som används i bruksanvisningen

Följ anvisningarna i bruksanvisningen för att inte utsätta dig själv och din omgivning för fara.

Allmän fara	<u>A</u>	Elrelaterad fara
Biologisk fara		Risk för skärskador
Fara orsakad av brännbara material	i	Avser information som är viktig, men inte riskrelaterad.
Risk för finger- eller handskador orsakade av mekaniska delar som oavsiktligt stängs.		

## Säkerhetsinstruktioner



Om inte dessa säkerhetsinstruktioner följs kan farliga situationer uppstå som kan resultera i dödsfall eller allvarliga personskador.

VARNING

- Följ säkerhetsföreskrifterna.
- Skakapparaten ska endast användas för avsedd användning. Felaktig användning kan leda till skador, kontaminering och personskador med dödlig utgång.
- Skakapparaten får endast användas av utbildad personal.
- Det är operatörens skyldighet att se till att lämplig personlig skyddsutrustning används. Användaren ska vara förtrogen med handboken "Laboratory Biosafety Manual" (utgiven av WHO) samt gällande nationella bestämmelser.



#### Skador från fel strömförsörjning.

Se till att skakapparaten är inkopplad endast i uttag som är ordentligt jordade.

VARNING



VARNING

#### Risker vid hantering av farliga ämnen.

Vid arbete med prover med korrosiva substanser (saltlösningar, syror, baser) måste tillbehören och skakapparaten rengöras grundligt.

- Skakapparaten är varken inert eller skyddad mot explosion. Använd aldrig skakapparaten i en miljö med fara för explosioner.
- Skaka inte giftiga eller radioaktiva ämnen eller patogena mikroorganismer utan lämpliga säkerhetsåtgärder.
- Vid skakning av farliga ämnen: Beakta Världshälsoorganisationen WHO:s handbok "Laboratory Biosafety Manual" och eventuella lokala bestämmelser. Vid skakning av mikrobiologiska prover från riskgrupp II (enligt Världshälsoorganisationen WHO:s handbok "Laboratory Biosafety Manual") måste aerosoltäta biologiska tätningar användas. Sök på internetsidan från Världshälsoorganisationen WHO (www.who.int) efter "Laboratory Biosafety Manual". För material i en högre riskgrupp måste extra säkerhetsåtgärder vidtas.
- Om toxiner eller patologiska substanser har kontaminerat skakapparaten eller dess delar måste lämpliga desinfektionsåtgärder och dekontamineringsmetoder vidtas ("Dekontaminering" på sidan 124, "Desinficering" på sidan 124).
- Om en farlig situation inträffar: Stäng av strömförsörjningen till skakapparaten och lämna området omedelbart.



VARNING

#### Hälsoskador från smittsamma ämnen.

Om vätskor eller andra material spills ut och hamnar under plattformen: Stäng omedelbart av skakapparaten, dra ur kontakten och ta bort plattformen ("Plattformar" på sidan 18). Städa upp spillet enligt normala laboratorierutiner. Använd lämplig personlig skyddsutrustning.



# Hälsorisker vid skakning av explosiva eller brännbara material eller ämnen.

Skaka inte explosiva eller brännbara material eller ämnen.

VARNING



#### Brännskador på heta ytor.

SE UPP



SE UPP

#### Skärskador från glasskärvor.

Om inte plattform och tillbehör, som kolvhållare, sitter fast ordentligt kan det leda till skadat glas om kärlen faller av skakapparaten.

I en temperaturkontrollerad skakapparat kan ytor såsom plattform och tillbehör under skakapparatens lock bli heta.

Vidrör inte heta ytor. Avvakta tills heta ytor har svalnat.

Se till att plattformen och tillbehören är korrekt monterade genom att använda rätt verktyg och skruvar. Se till att kolvhållarna klarar belastningen av kärlet och provet vid vald hastighet. Kolvhållare med din individuellt konfigurerade belastning kan ha lägre hastighetskapacitet än hållfasthetsbegränsningen hos skakapparaten. Se "3. 4. Lastning och normal användning" på sidan 115. Var uppmärksam på ovanliga ljud. De kan indikera att plattform eller tillbehör inte sitter fast.



SF IIPP

## Skador på skakapparaten eller felfunktion till följd av skadad pekskärm.

- Använd inte skakapparaten om pekskärmen är skadad.
- Stäng av skakapparaten. Koppla ur stickkontakten. Få pekskärmen utbytt av en auktoriserad service tekniker.



**Biologisk skada orsakad av trasiga eller läckande kärl.** Felaktigt monterade tillbehör kan leda till att prover spills ut.

SE UPP

- Se till att tillbehören är korrekt monterade genom att använda rätt verktyg och skruvar.
- Se till att tillbehören passar på plattformen.
- Använd alltid kärl med tillbehör av passande storlek.
- Kärlen måste vara intakta och monterade på rätt sätt.



Säkerheten kan försämras genom felaktig lastning och skadade tillbehör.

- SE UPP
- Se alltid till att lasten (tillbehör och prover) är fördelad så jämnt som möjligt, särskilt när du använder en dual stack-plattform.
- Använd inte tillbehör som uppvisar tecken på korrosion eller sprickor. Kontakta kundtjänst för mer information.
- Använd endast skakapparater som har lastats korrekt.
- Overbelasta aldrig skakapparaten.
- Kontrollera att tillbehören är korrekt monterade innan du använder skakapparaten. Följ instruktionerna i avsnitt "Tillbehör" på sidan 88.



SE UPP

Fysisk skada orsakad av att de grundläggande användningsrutinerna åsidosätts.

- Använd aldrig skakapparaten utan en korrekt monterad plattform.
- Använd aldrig skakapparaten om delar av utsidan är skadade eller saknas.
- Använd aldrig en skakapparat med skadad lockfjäder. En defekt lockfjäder kan inte hålla locket i fullt öppet läge på ett säkert sätt.
- Flytta inte skakapparaten medan den är igång.
- Luta dig inte mot skakapparaten.
- Fyll eller töm aldrig skakapparaten förrän den har stannat . helt och detta har bekräftats på pekskärmen.
- Placera inte något på skakapparaten medan den är igång.
- Vidrör inte plattformen eller några tillbehör på skakapparaten när den är igång.
- Skakapparatens kåpa får inte öppnas av operatören.



OBS

#### Skyddet kan försämras av inkompatibla tillbehör.

Använd endast tillbehör för skakapparaten, som har godkänts av Thermo Fisher Scientific. Aktuella listor finns på www. thermofisher.com.



OBS

#### Så här stänger du av skakapparaten:

Tryck på knappen STOPP. Stäng av skakapparaten med huvudströmbrytaren. Dra ur stickproppen. Om ett nödläge uppstår, frånkoppla strömförsörjningen.

## 1. Specifikationer

## 1.1. Tekniska data

Thermo Scientific Solaris 2000 skakapparat					
Hastighetsområde	15 – 525 v/min				
Körtid	99 h 59 min (1 min inter	vall) eller kontinuerligt läge			
Bullemivå vid max. hastighet	50 dB (A) (1 m framför e	nheten vid 1,6 m höjd)			
Maximal belastning (inkl. plattform, tillbehör och prover)	25 kg (55 lbs)				
Elektrisk anslutning	100–240 V ±10 %, 50 /	' 60 Hz			
Effektförbrukning	60 W				
Krav på omgivningen					
För förvaring och transport	Temperatur: -10 °C till 5 Fuktighet: 15 % till 85 %	5 °C			
För användning	Använd inomhus Altituder upp till 3 000 meter över havet Temperatur: 5 °C till 40 °C Max. relativ luftfuktighet 80 % upp till 31 °C; linjärt avtagande ner till 50 % relativ luftfuktighet vid 40 °C				
Nedsmutsningsgrad	2				
Överspänningskategori	II				
IP	20				
Gränssnitt	USB	Ethernet			
	2x USB-A 2.0	RJ45			
Dimensioner	Längd	Bredd	Höjd		
	47 cm (18,5 in)	37 cm (14,5 in)	15 cm (5,5 in)		
Vikt	20,9 kg (46,0 lbs)				

Tabell 1: Tekniska data Solaris 2000

Thermo	Scientific	Solaris	4000	Skakappara	at
--------	------------	---------	------	------------	----

Hastighetsområde	15 – 525 v/min			
Körtid	99 h 59 min (1 min intervall) eller kontinuerligt läge			
Bullemivå vid max. hastighet	54 dB (A) (1 m framför e	nheten vid 1,6 m höjd)		
Maximal belastning (inkl. plattform, tillbehör och prover)	43 kg (95 lbs)			
Elektrisk anslutning	100–240 V ±10 %, 50 /	60 Hz		
Effektförbrukning	80 W			
Krav på omgivningen				
För förvaring och transport	Temperatur: -10 °C till 5 Fuktighet: 15 % till 85 %	5 °C		
För användning	Använd inomhus Altituder upp till 3 000 meter över havet Temperatur: 5 °C till 40 °C Max. relativ luftfuktighet 80 % upp till 31 °C; linjärt avtagande ner till 50 % relativ luftfuktighet vid 40 °C			
Nedsmutsningsgrad	2			
Överspänningskategori	11			
P	20			
Gränssnitt	USB	Ethernet		
	2x USB-A 2.0	RJ45		
Dimensioner	Längd	Bredd	Höjd	
	65 cm (25,6 in)	58 cm (22,8 in)	18 cm (7 in)	
Vikt	75,1 kg (165,5 lbs)			

Tabell 2: Tekniska data Solaris 4000

## Thermo Scientific Solaris 2000 I skakapparat

Hastighetsområde	15 – 525 v/min		
Temperaturområde	30-60 °C		
Temperaturstabilitet i kolv vid 37 °C	Omgivningstemperatur 2 ± 0,1 °C	3 °C; i 1 timme med stabi	kammartemperatur
Homogen temperatur i kolv vid 37 °C	Omgivningstemperatur 2 ± 0,5 °C	3 °C; i 1 timme med stabi	kammartemperatur
Körtid	99 h 59 min (1 min interv	vall) eller kontinuerligt läge	
Bullernivå vid max. hastighet	52 dB (A) (1 m framför er	nheten vid 1,6 m höjd)	
Maximal belastning (inkl. plattform, tillbehör och prover)	16 kg (35 lbs), inklusive 11 x 14-plattfor	m på 2,3 kg (5,1 lbs)	
Elektrisk anslutning	100–120, 200–240 V ±	10 %, 50 / 60 Hz	
Effektförbrukning	900 W		
Krav på omgivningen			
För förvaring och transport	Temperatur: -10 °C till 55 °C Fuktighet: 15 % till 85 %		
För användning	Använd inomhus Altituder upp till 3 000 meter över havet Temperatur: 5 °C till 40 °C Max. relativ luftfuktighet 80 % upp till 31 °C; linjärt avtagande ner till 50 % relativ luftfuktighet vid 40 °C		
Nedsmutsningsgrad	2		
Överspänningskategori	1		
IP	20		
Gränssnitt	USB	Ethernet	
	2x USB-A 2.0	RJ45	
Dimensioner	Längd	Bredd	Höjd
	70 cm (27,6 in)	36 cm (14,2 in)	46 cm (18,1 in)
Med öppet lock			79 cm (30,8 in)
Vikt	44,9 kg (99 lbs)		

Tabell 3: Tekniska data Solaris 2000 I

## Thermo Scientific Solaris 4000 I skakapparat

Hastighetsområde	15 – 525 v/min			
Temperaturområde	30–60 °C			
Temperaturstabilitet i kolv vid 37 °C	Omgivningstemperatur 2 ± 0,1 °C	Omgivningstemperatur 23 °C; i 1 timme med stabil kammartemperatur $\pm$ 0,1 °C		
Homogen temperatur i kolv vid 37 °C	Omgivningstemperatur 2 ± 0,5 °C	3 °C; i 1 timme med stabi	l kammartemperatur	
Körtid	99 h 59 min (1 min inter	vall) eller kontinuerligt läge		
Bullemivå vid max. hastighet	52 dB (A) (1 m framför e	nheten vid 1,6 m höjd)		
Maximal belastning (inkl. plattform, tillbehör och prover)	32 kg (70,5 lbs), inklusive 18 x 18-plattfor	m på 4,3 kg (9,5 lbs)		
Elektrisk anslutning	100–120, 200–240 V ±	10 %, 50 / 60 Hz		
Effektförbrukning	900 W			
Krav på omgivningen				
För förvaring och transport	Temperatur: -10 °C till 55 °C Fuktighet: 15 % till 85 %			
För användning	Använd inomhus Altituder upp till 3 000 meter över havet Temperatur: 5 °C till 40 °C Max. relativ luftfuktighet 80 % upp till 31 °C; linjärt avtagande ner till 50 % relativ luftfuktighet vid 40 °C			
Nedsmutsningsgrad	2			
Överspänningskategori	11			
IP	20			
Gränssnitt	USB	Ethernet		
	2x USB-A 2.0	RJ45		
Dimensioner	Längd	Bredd	Höjd	
	77 cm (30,3 in)	57 cm (22,4 in)	55 cm (21,7 in)	
Med öppet lock			94 cm (36,7 in)	
Vikt	68,1 kg (150,1 lbs)			

Tabell 4: Tekniska data Solaris 4000 I

## Thermo Scientific Solaris 2000 R skakapparat

Hastighetsområde	15 – 525 v/min			
Temperaturområde	5–60 °C	5–60 °C		
Temperaturstabilitet i kolv vid 37 °C	Omgivningstemperatur 23 °C; i 1 timme med stabil kammartemperatur $\pm$ 0,1 °C			
Homogen temperatur i kolv vid 37 °C	Omgivningstemperatur 2 ± 0,5 °C	3 °C; i 1 timme med stabi	kammartemperatur	
Körtid	99 h 59 min (1 min interv	vall) eller kontinuerligt läge		
Bullernivå vid max. hastighet	56 dB (A) (1 m framför er	nheten vid 1,6 m höjd)		
Maximal belastning (inkl. plattform, tillbehör och prover)	16 kg (35 lbs), inklusive 11 x 14-plattfor	m på 2,3 kg (5,1 lbs)		
Elektrisk anslutning	100–240 V ±10 %, 50 /	60 Hz		
Effektförbrukning	350 W			
Krav på omgivningen				
För förvaring och transport	Temperatur: -10 °C till 55 °C Fuktighet: 15 % till 85 %			
För användning	Använd inomhus Altituder upp till 3 000 meter över havet Temperatur: 5 °C till 40 °C Max. relativ luftfuktighet 80 % upp till 31 °C; linjärt avtagande ner till 50 % relativ luftfuktighet vid 40 °C			
Nedsmutsningsgrad	2			
Överspänningskategori	l			
IP	20			
Gränssnitt	USB	Ethernet		
	2x USB-A 2.0	RJ45		
Dimensioner	Längd	Bredd	Höjd	
	70 cm (27,6 in)	36 cm (14,2 in)	46 cm (18,1 in)	
Med öppet lock			79 cm (30,8 in)	
Vikt	47,2 kg (104,1 lbs)			

Tabell 5: Tekniska data Solaris 2000 R

## Thermo Scientific Solaris 4000 R skakapparat

Hastighetsområde	15 – 525 v/min			
Temperaturområde	4–60 °C	4–60 °C		
Temperaturstabilitet i kolv vid 37 °C	Omgivningstemperatur 23 °C; i 1 timme med stabil kammartemperatur $\pm$ 0,1 °c			
Homogen temperatur i kolv vid 37 °C	Omgivningstemperatur 23 ± 0,5 °C	3 °C; i 1 timme med stabil	kammartemperatur	
Körtid	99 h 59 min (1 min interva	all) eller kontinuerligt läge		
Bullernivå vid max. hastighet	56 dB (A) (1 m framför en	heten vid 1,6 m höjd)		
Maximal belastning (inkl. plattform, tillbehör och prover)	32 kg (70,5 lbs), inklusive 18 x 18-plattform	n på 4,3 kg (9,5 lbs)		
Elektrisk anslutning	100–240 V ±10 %, 50 / 6	60 Hz		
Effektförbrukning	750 W			
Krav på omgivningen				
För förvaring och transport	Temperatur: -10 °C till 55 °C Fuktighet: 15 % till 85 %			
För användning	Använd inomhus Altituder upp till 3 000 meter över havet Temperatur: 5 °C till 40 °C Max. relativ luftfuktighet 80 % upp till 31 °C; linjärt avtagande ner till 50 % relativ luftfuktighet vid 40 °C			
Nedsmutsningsgrad	2			
Överspänningskategori	1			
P	20			
Gränssnitt	USB	Ethernet		
	2x USB-A 2.0	RJ45		
Dimensioner	Längd	Bredd	Höjd	
	77 cm (30,3 in)	57 cm (22,4 in)	55 cm (21,7 in)	
Med öppet lock			94 cm (36,7 in)	
Vikt	74,9 kg (165,1 lbs)			

Tabell 6: Tekniska data Solaris 4000 R

## 1.2. Tillbehör



Skyddet kan försämras av inkompatibla tillbehör. Använd endast tillbehör för skakapparaten, som har godkänts av Thermo Fisher Scientific.

Aktuella listor finns på <u>www.thermofisher.com</u>.

## 1.2.1. Plattformar

Plattform		
Solaris 2000	Vikt	
Thermo Scientific Solaris 12 x 14 universalplattform	2,7 kg (5,9 lbs)	SK1214
Thermo Scientific Solaris 12 x 14 Dual Stack universalplattform	6,4 kg (14,1 lbs)	SK1214D
Thermo Scientific Solaris 12 x 14 Dual stack universalplattform, uppgraderingssats		SK1214DK
Thermo Scientific Solaris 18 x 18 universalplattform	4,3 kg (9,4 lbs)	SK1818
Thermo Scientific Solaris 18 x 18 Dual Stack universalplattform	9,7 kg (21,3 lbs)	SK1818D
Thermo Scientific Solaris 18 x 18 Dual stack universalplattform, uppgraderingssats		SK1818DK
Thermo Scientific Solaris 18 x 24 universalplattform	6,0 kg (13,2 lbs)	SK1824
Solaris 4000		
Thermo Scientific Solaris 18 x 30 universalplattform	7,5 kg (16,5 lbs)	SK1830
Thermo Scientific Solaris 18 x 30 Dual Stack universalplattform	16,3 kg (35,9 lbs)	SK1830D
Thermo Scientific Solaris 18 x 30 Dual stack universalplattform, uppgraderingssats		SK1830DK
Thermo Scientific Solaris 36 x 24 universalplattform	12,0 kg (26,4 lbs)	SK3624
Solaris 2000 I / 2000 R		
Thermo Scientific Solaris 11 x 14 universalplattform	2,3 kg (5,1 lbs)	SK1114
Solaris 4000 I / 4000 R		
Thermo Scientific Solaris 18 x 18 universalplattform	4,3 kg (9,5 lbs)	SK1818
Reservdelar och tillbehör		
Klämma, Reservdelssats (Skruvar)		SK1001
Liten reservdelssats för plattform (plattformsskruvar, verktyg)		SK0100
Stor reservdelssats för plattform (plattformsskruvar, verktyg)		SK0101
Skruvmejsel för tillbehör		75004131

Tabell 7: Tillgängliga plattformar

## 1.2.2. Kolvhållare



## För Solaris 2000-plattformar

	Artikel- nummer	Universal			Dual	Stack
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
10 ml Erlenmeyer	30150BI	72	113	157	140	226
25 ml Erlenmeyer	30151	42	64	80	80	124
50 ml Erlenmeyer	30152BI	42	64	80	80	124
125 ml Erlenmeyer	30153	15	32	40	30	62
250 ml Erlenmeyer	30154BI	9	16	24	16	32
300 ml Erlenmeyer	30155	9	16	20	16	32
500 ml Erlenmeyer	30156BI	9	16	20	16	32
1 l Erlenmeyer	30157BI	4	9	11	8	16
2 l Erlenmeyer	30158	3	5	6	-	-
4 l Erlenmeyer	30159	1	4	4	-	-
5 l Erlenmeyer	30159B	1	2	4	-	-
6 l Erlenmeyer	30160	1	2	2	-	-
2800 ml Fernbach kolv	30162	1	4	4	-	-
Låg odlingskolv 2,5 l	30161	1	2	2	_	_

Tabell 8: Tillgängliga kolvhållare för Solaris 2000-plattform

## För Solaris 4000-plattformar

	Artikel- nummer	Universal		Dual Stack
		18 x 30	36 x 24	18 x 30
10 ml Erlenmeyer	30150BI	203	187	402
25 ml Erlenmeyer	30151	112	187	220
50 ml Erlenmeyer	30152BI	112	187	220
125 ml Erlenmeyer	30153	46	83	92
250 ml Erlenmeyer	30154BI	34	40	68
300 ml Erlenmeyer	30155	28	40	56
500 ml Erlenmeyer	30156BI	28	40	56
1   Erlenmeyer	30157BI	14	20	28
2 l Erlenmeyer	30158	8	11	16
4 I Erlenmeyer	30159	6	8	-
5 l Erlenmeyer	30159B	5	8	-
6 l Erlenmeyer	30160	3	6	-
2800 ml Fernbach kolv	30162	6	8	12
Låg odlingskolv 2,5 l	30161	3	6	6

Tabell 9: Tillgängliga kolvhållare för Solaris 4000-plattform

## För Solaris 2000 I / 2000 R-plattformar

	Artikel- nummer	Universal
		11 x 14
10 ml Erlenmeyer	30150BI	59
25 ml Erlenmeyer	30151	35
50 ml Erlenmeyer	30152BI	35
125 ml Erlenmeyer	30153	15
250 ml Erlenmeyer	30154BI	7
300 ml Erlenmeyer	30155	7
500 ml Erlenmeyer	30156BI	7
1   Erlenmeyer	30157BI	4
2 l Erlenmeyer	30158	2
4 l Erlenmeyer	30159	-
5 l Erlenmeyer	30159B	-
6 l Erlenmeyer	30160	-
2800 ml Fernbach kolv	30162	1
Låg odlingskolv 2,5 l	30161	1

Tabell 10: Tillgängliga kolvhållare för Solaris 2000 I/2000 R-plattform

För	Solaris	4000	17	4000	<b>R-plattformar</b>

	Artikel- nummer	Universal
		18 x 18
10 ml Erlenmeyer	30150BI	113
25 ml Erlenmeyer	30151	64
50 ml Erlenmeyer	30152BI	64
125 ml Erlenmeyer	30153	32
250 ml Erlenmeyer	30154BI	16
300 ml Erlenmeyer	30155	16
500 ml Erlenmeyer	30156BI	16
1   Erlenmeyer	30157BI	9
2 l Erlenmeyer	30158	6
4 I Erlenmeyer	30159	-
5 l Erlenmeyer	30159B	-
6 l Erlenmeyer	30160	-
2800 ml Fernbach kolv	30162	4
Låg odlingskolv 2,5 l	30161	1

Tabell 11: Tillgängliga kolvhållare för Solaris 4000 I/4000 R-plattform

## 1. 2. 3. Hållare för mikrotiterplatta/deepwellplatta



## För Solaris 2000-plattformar

	Artikel- nummer		Universal		Dual	Stack
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Mikrotiterplatta / Deepwellplatta	30175	7	12	16	14	24

Tabell 12: Kapacitet för hållare för mikrotiterplatta/deepwellplatta för Solaris 2000-plattform

## För Solaris 2000 I / 2000 R-plattformar

	Artikel- nummer	Universal
		11 x 14
Mikrotiterplatta / Deepwellplatta	30175	6

Tabell 13: Kapacitet för hållare för mikrotiterplatta/deepwellplatta för Solaris 2000 l/2000R-plattform

#### För Solaris 4000-plattformar

	Artikel- nummer	Universal		Dual Stack
		18 x 30	36 x 24	18 x 30
Mikrotiterplatta / Deepwellplatta	30175	23	36	42

Tabell 14: Kapacitet för hållare för mikrotiterplatta/deepwellplatta för Solaris 4000-plattform

#### För Solaris 4000 I / 4000 R-plattformar

	Artikel- nummer	Universal
		18 x 18
Mikrotiterplatta / Deepwellplatta	30175	10

Tabell 15: Kapacitet för hållare för mikrotiterplatta/deepwellplatta för Solaris 4000 I/4000R-plattform

## 1. 2. 4. Provrörsställ

#### För Solaris 2000-plattformar



	Artikel- nummer	Universal		Dual Stack		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Halv storlek						
10 – 13 mm, röd, 6 x 6 hål	30181	8	12	15	14	22
14 – 16 mm, orange, 6 x 6 hål	30183	5	9	11	9	16
17 – 20 mm, vit, 4 x 5 hål	30185	7	11	14	12	20
21 – 25 mm, blå, 4 x 4 hål	30187	6	9	11	10	16
26 – 30 mm, grön, 3 x 3 hål	30189	6	9	12	11	17
Mikrocentrifug, 1,5 ml, blå, 4 x 6 hål	30191	6	10	13	11	19
Full storlek						
10 – 13 mm, 6 x 12 hål	30180BI	З	7	9	6	12
14 – 16 mm, 6 x 12 hål	30182	3	4	6	6	8
17 – 20 mm, 4 x 10 hål	30184	3	5	7	6	8
21 – 25 mm, 4 x 10 hål	30186	2	3	5	4	6
26 – 30 mm, 3 x 8 hål	30188	3	4	5	5	8
1,5 ml mikrocentrifug, 8 x 12 hål	30190	3	4	7	6	8

Tabell 16: Tillgängliga provrörsställ för Solaris 2000 plattformar

## För Solaris 4000-plattformar

	Artikel- nummer	Universal		Dual Stack
		18 x 30	36 x 24	18 x 30
Halv storlek				
10 – 13 mm, röd, 6 x 6 hål	30181	21	32	40
14 – 16 mm, orange, 6 x 6 hål	30183	13	20	24
17 – 20 mm, vit, 4 x 5 hål	30185	18	20	34
21 – 25 mm, blå, 4 x 4 hål	30187	14	22	26
26 – 30 mm, grön, 3 x 3 hål	30189	15	24	30
1,5 ml mikrocentrifug, blå, 4 x 6 hål	30191	18	24	34
Full storlek				
10 – 13 mm, 6 x 12 hål	30180BI	13	20	18
14 – 16 mm, 6 x 12 hål	30182	9	12	16
17 – 20 mm, 4 x 10 hål	30184	9	15	18
21 – 25 mm, 4 x 10 hål	30186	6	9	12
26 – 30 mm, 3 x 8 hål	30188	7	10	13
1,5 ml mikrocentrifug, 8 x 12 hål	30190	9	12	17

Tabell 17: Tillgängliga provrörsställ för Solaris 4000 plattformar

## För Solaris 2000 I / 2000 R-plattformar

	Artikel- nummer	Universal
		11 x 14
Halv storlek		
10 – 13 mm, röd, 6 x 6 hål	30181	6
14 – 16 mm, orange, 6 x 6 hål	30183	2
17 – 20 mm, vit, 4 x 5 hål	30185	4
21 – 25 mm, blå, 4 x 4 hål	30187	2
26 – 30 mm, grön, 3 x 3 hål	30189	4
Mikrocentrifug, 1,5 ml, blå, 4 x 6 hål	30191	3
Full storlek		
10 – 13 mm, 6 x 12 hål	30180BI	3
14 – 16 mm, 6 x 12 hål	30182	2
17 – 20 mm, 4 x 10 hål	30184	2
21 – 25 mm, 4 x 10 hål	30186	1
26 – 30 mm, 3 x 8 hål	30188	2
1,5 ml mikrocentrifug, 8 x 12 hål	30190	2

Tabell 18: Tillgängliga provrörsställ för Solaris 2000 I / 2000 R plattformar

## För Solaris 4000 I / 4000 R-plattformar

	Artikel- nummer	Universal
		18 x 18
Halv storlek		
10 – 13 mm, röd, 6 x 6 hål	30181	12
14 – 16 mm, orange, 6 x 6 hål	30183	8
17 – 20 mm, vit, 4 x 5 hål	30185	10
21 – 25 mm, blå, 4 x 4 hål	30187	8
26 – 30 mm, grön, 3 x 3 hål	30189	9
Mikrocentrifug, 1,5 ml, blå, 4 x 6 hål	30191	8
Full storlek		
10 – 13 mm, 6 x 12 hål	30180BI	7
14 – 16 mm, 6 x 12 hål	30182	4
17 – 20 mm, 4 x 10 hål	30184	5
21 – 25 mm, 4 x 10 hål	30186	3
26 – 30 mm, 3 x 8 hål	30188	4
1,5 ml mikrocentrifug, 8 x 12 hål	30190	5

Tabell 19: Tillgängliga provrörsställ för Solaris 4000 I / 4000 R plattformar

## 1. 2. 5. Provrörsställhållare med justerbar vinkel

### För Solaris 2000-plattformar



	Artikel- nummer	Universal		Dual Stack		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
10–13 mm, 72 platser	236090	3	4	6	6	8
16–20 mm, 40 platser	236091	3	4	6	5	8
21–25 mm, 40 platser	236092	3	3	4	5	6
26–30 mm, 24 platser	236093	3	3	4	5	6

Tabell 20: Tillgängliga provrörsställhållare med justerbar vinkel för Solaris 2000-plattform

## För Solaris 4000-plattformar

	Artikel- nummer	Universal		Dual Stack
		18 x 30	36 x 24	18 x 30
10–13 mm, 72 platser	236090	7	14	14
16–20 mm, 40 platser	236091	7	14	14
21–25 mm, 40 platser	236092	5	9	10
26–30 mm, 24 platser	236093	6	9	12

Tabell 21: Tillgängliga provrörsställhållare med justerbar vinkel för Solaris 4000-plattform

## För Solaris 2000 I / 2000 R-plattformar

	Artikel- nummer	Universal
		11 x 14
10–13 mm, 72 platser	236090	2
16–20 mm, 40 platser	236091	2
21–25 mm, 40 platser	236092	1
26–30 mm, 24 platser	236093	1

Tabell 22: Tillgängliga provrörsställhållare med justerbar vinkel för Solaris 2000 I/2000 R-plattform

## För Solaris 4000 I / 4000 R-plattformar

	Artikel- nummer	Universal
		18 x 18
10–13 mm, 72 platser	236090	4
16–20 mm, 40 platser	236091	3
21–25 mm, 40 platser	236092	3
26–30 mm, 24 platser	236093	3

Tabell 23: Tillgängliga provrörsställhållare med justerbar vinkel för Solaris 4000 I/4000 R-plattform

## 1. 2. 6. Justerbara kärlhållare



#### För Solaris 2000-plattformar

	Artikel- nummer	Universal		Dual	Stack	
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Kärlhållare justerbar i höjd, 11" x 14"	75004104	-	-	-	-	-
Kärlhållare justerbar i två höjder, 12" x 14"	75004102	1	1	1	-	2
Kärlhållare justerbar i två höjder, 18" x 18"	75004103	-	1	1	-	-
Justerbar kärlhållare	75004101	4	4	6	6	8

Tabell 24: Tillgängliga justerbara kärlhållare för Solaris 2000-plattform

## För Solaris 4000-plattformar

	Artikel- nummer	Unive	ersal	Dual Stack
		18 x 30	36 x 24	18 x 30
Kärlhållare justerbar i höjd, 11" x 14"	75004104	_	_	_
Kärlhållare justerbar i två höjder, 12" x 14"	75004102	2	2	4
Kärlhållare justerbar i två höjder, 18" x 18"	75004103	1	2	2
Justerbar kärlhållare	75004101	8	12	16

Tabell 25: Justerbara kärlhållare för Solaris 4000-plattform

## För Solaris 2000 I / 2000 R-plattformar

	Artikel- nummer	Universal
		11 x 14
Kärlhållare justerbar i höjd, 11" x 14"	75004104	1
Kärlhållare justerbar i två höjder, 12" x 14"	75004102	-
Kärlhållare justerbar i två höjder, 18" x 18"	75004103	-
Justerbar kärlhållare	75004101	2

Tabell 26: Justerbara kärlhållare för Solaris 2000 I/2000 R-plattform

## För Solaris 4000 I / 4000 R-plattformar

	Artikel- nummer	Universal
		18 x 18
Kärlhållare justerbar i höjd, 11" x 14"	75004104	-
Kärlhållare justerbar i två höjder, 12" x 14"	75004102	1
Kärlhållare justerbar i två höjder, 18" x 18"	75004103	1
Justerbar kärlhållare	75004101	4

Tabell 27: Justerbara kärlhållare för Solaris 4000 I/4000 R-plattform

## 1. 2. 7. Kvadratiska mediahållare

## För Solaris 2000-plattformar



	Artikel- nummer	Universal		Dual Stack		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
125 ml kvadratisk kolvhållare	75004106	12	16	24	21	32
250 ml kvadratisk kolvhållare	75004107	9	16	22	18	32
500 ml kvadratisk kolvhållare	75004108	6	9	12	11	18
1 000 ml kvadratisk kolvhållare	75004109	4	9	12	8	18

Tabell 28: Tillgängliga kvadratiska mediahållare för Solaris 2000-plattform

## För Solaris 4000-plattformar

	Artikel- nummer	Univ	ersal	Dual Stack	
		18 x 30	36 x 24	18 x 30	
125 ml kvadratisk kolvhållare	75004106	28	54	56	
250 ml kvadratisk kolvhållare	75004107	28	52	56	
500 ml kvadratisk kolvhållare	75004108	15	24	30	
1 000 ml kvadratisk kolvhållare	75004109	15	24	28	

Tabell 29: Tillgängliga kvadratiska mediahållare för Solaris 4000-plattform

## För Solaris 2000 I / 2000 R-plattformar

	Artikel- nummer	Universal
		11 x 14
125 ml kvadratisk kolvhållare	75004106	9
250 ml kvadratisk kolvhållare	75004107	7
500 ml kvadratisk kolvhållare	75004108	3
1 000 ml kvadratisk kolvhållare	75004109	2

Tabell 30: Tillgängliga kvadratiska mediahållare för Solaris 2000 I/2000 R-plattform

## För Solaris 4000 I / 4000 R-plattformar

	Artikel- nummer	Universal
		18 x 18
125 ml kvadratisk kolvhållare	75004106	16
250 ml kvadratisk kolvhållare	75004107	16
500 ml kvadratisk kolvhållare	75004108	9
1 000 ml kvadratisk kolvhållare	75004109	9

Tabell 31: Tillgängliga kvadratiska mediahållare för Solaris 4000 I/4000 R-plattform

## 1. 2. 8. Nalgene bägarställ

### För Solaris 2000-plattformar



Tabell 32: Tillgängliga Nalgene bägarställ för Solaris 2000-plattform



## För Solaris 4000-plattformar

	Artikel- nummer	Universal		Dual Stack
		18 x 30	36 x 24	18 x 30
Storlek 11 x 14/12 x 14			Ställ/platser	
Nalgene 30 ml bägare, 36 platser	75004116	-	-	-
Nalgene 50 ml bägare, 28 platser	75004129	-	-	-
Nalgene 100 ml bägare, 15 platser	75004118	-	-	-
Nalgene 250 ml bägare, 9 platser	75004119	-	-	-
Nalgene 400 ml bägare, 8 platser	75004120	-	-	-
Nalgene 600 ml bägare, 5 platser	75004121	-	-	-
Storlek 9 x 18			Ställ/platser	
Nalgene 30 ml bägare, 32 platser	75004110	3/96	4 / 128	6 / 192
Nalgene 50 ml bägare, 24 platser	75004128	3/72	4 / 96	6 / 144
Nalgene 100 ml bägare, 15 platser	75004112	3 / 45	4 / 60	6/90
Nalgene 250 ml bägare, 8 platser	75004113	3/24	4 / 32	6/48
Nalgene 400 ml bägare, 6 platser	75004114	3/18	4 / 24	6/36
Nalgene 600 ml bägare, 4 platser	75004115	3/12	4 / 16	6 / 24

Tabell 33: Tillgängliga Nalgene bägarställ för Solaris 4000-plattform

## För Solaris 2000 I / 2000 R-plattformar

	Artikel- nummer	Universal
Storlek 11 x 14/12 x 14		Ställ/platser
Nalgene 30 ml bägare, 36 platser	75004116	1/36
Nalgene 50 ml bägare, 28 platser	75004129	1 / 28
Nalgene 100 ml bägare, 16 platser	75004118	1 / 16
Nalgene 250 ml bägare, 9 platser	75004119	1/9
Nalgene 400 ml bägare, 8 platser	75004120	1/8
Nalgene 600 ml bägare, 5 platser	75004121	1/5
Storlek 9 x 18		Ställ/platser
Nalgene 30 ml bägare, 32 platser	75004110	-
Nalgene 50 ml bägare, 24 platser	75004128	-
Nalgene 100 ml bägare, 15 platser	75004112	-
Nalgene 250 ml bägare, 8 platser	75004113	-
Nalgene 400 ml bägare, 6 platser	75004114	_
Nalgene 600 ml bägare, 4 platser	75004115	-

Tabell 34: Tillgängliga Nalgene bägarställ för Solaris 2000 I/2000 R-plattform
Tillbehör

### För Solaris 4000 I / 4000 R-plattformar

	Artikel- nummer	Universal
		18 x 18
Storlek 11 x 14/12 x 14		Ställ/platser
Nalgene 30 ml bägare, 36 platser	75004116	-
Nalgene 50 ml bägare, 28 platser	75004129	-
Nalgene 100 ml bägare, 16 platser	75004118	-
Nalgene 250 ml bägare, 9 platser	75004119	-
Nalgene 400 ml bägare, 8 platser	75004120	-
Nalgene 600 ml bägare, 5 platser	75004121	-
Storlek 9 x 18		Ställ/platser
Nalgene 30 ml bägare, 32 platser	75004110	2/64
Nalgene 50 ml bägare, 24 platser	75004128	2/48
Nalgene 100 ml bägare, 15 platser	75004112	2/30
Nalgene 250 ml bägare, 8 platser	75004113	2/16
Nalgene 400 ml bägare, 6 platser	75004114	2/12
Nalgene 600 ml bägare, 4 platser	75004115	2/8

Tabell 35: Tillgängliga Nalgene bägarställ för Solaris 4000 I/4000 R-plattform

### 1. 2. 9. Separertratthållare

### För Solaris 2000-plattformar



	Artikel- nummer	Universal		Dual	Stack	
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Separertratthållare, 250 ml till 2 000 ml	75004125	2	З	4	2	5
Vertikalt monterad separertratthållare		4	4	6	-	-

 Tabell 36:
 Tillgängliga separertratthållare för Solaris 2000-plattform

### För Solaris 4000-plattformar

	Artikel- nummer	Unive	ersal	Dual Stack
		18 x 30	36 x 24	18 x 30
Separertratthållare, 250 ml till 2000 ml	75004125	5	7	9
Vertikalt monterad separertratthållare		8	12	-

 Tabell 37: Tillgängliga separertratthållare för Solaris 4000-plattform

### För Solaris 2000 I / 2000 R-plattformar

	Artikelnummer	Universal
Separertratthållare, 250 ml till 2000 ml	75004125	-
Vertikalt monterad separertratthållare		-

Tabell 38: Tillgängliga separertratthållare för Solaris 2000 I/2000 R-plattform

#### För Solaris 4000 I / 4000 R-plattformar

	Artikelnummer	Universal
		18 x 18
Separertratthållare, 250 ml till 2000 ml	75004125	-
Vertikalt monterad separertratthållare		-

 Tabell 39: Tillgängliga separertratthållare för Solaris 4000 I/4000 R-plattform

### 1. 2. 10. Universalbricka

### För Solaris 2000-plattformar



	Artikelnummer	Universal		rtikelnummer Universal Dual Stack			Stack
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18	
11 x 14/12 x 14 universalbricka	75004123	1	-	-	2	-	
9 x 18 universalbricka	75004122	-	2	3	-	4	

Tabell 40: Tillgängliga universalbrickor för Solaris 2000-plattform

### För Solaris 4000-plattformar

	Artikelnummer Universal		ersal	Dual Stack
		18 x 30	36 x 24	18 x 30
11 x 14/12 x 14 universalbricka	75004123	-	-	-
9 x 18 universalbricka	75004122	3	6	6

 Tabell 41: Tillgängliga universalbrickor för Solaris 4000-plattform

### För Solaris 2000 I / 2000 R-plattformar

	Artikelnummer	Universal
		11 x 14
11 x 14/12 x 14 universalbricka	75004123	1
9 x 18 universalbricka	75004122	-

 Tabell 42: Tillgängliga universalbrickor för Solaris 2000 I/2000 R-plattform

### För Solaris 4000 I / 4000 R-plattformar

	Artikelnummer	Universal
		18 x 18
11 x 14/12 x 14 universalbricka	75004123	-
9 x 18 universalbricka	75004122	2

Tabell 43: Tillgängliga universalbrickor för Solaris 4000 I/4000 R-plattform

### Tillbehör för universalbrickor

	Artikelnummer	Universal
O-ringsset (universalbricka)	75004132	-

Tabell 44: Tillgängliga tillbehör för universalbrickor

### 1. 2. 11. Klistermattor

Allmänn information för val av klistermatta:

 Mattstorleken 170 x 280 mm är optimerad för plattformarna 11 x 14 och 12 x 14.



Alla mattor kan skäras till i önskad storlek.

#### För Solaris 2000-plattformar

	Artikel- nummer	Universal		tikel- Universal I mmer u I		Dual S univers plattfo	tack sal- rm
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18	
200 x 200 mm matta med hög vidhäftning, 2-pack	75004126	1	4	6	2	8	
280 x 170 mm matta med hög vidhäftning, 2-pack	75004127	2	2	4	4	4	
200 x 200 mm matta med låg vidhäftning, 2-pack	75004111	1	4	6	2	8	
280 x 170 mm matta med låg vidhäftning, 2-pack	75004117	2	2	4	4	4	

Mattorna säljs i förpackningar om 2. Antal mattor för en plattform avser inte denna förpackning, utan en enstaka matta.

 Tabell 45: Tillgängliga klistermattor för Solaris 2000-plattform

### För Solaris 4000-plattformar

	Artikel- nummer	Universal		Dual Stack
		18 x 30	36 x 24	18 x 30
200 x 200 mm matta med hög vidhäftning, 2-pack	75004126	6	12	12
280 x 170 mm matta med hög vidhäftning, 2-pack	75004127	6	9	12
200 x 200 mm matta med låg vidhäftning, 2-pack	75004111	6	12	12
280 x 170 mm matta med låg vidhäftning, 2-pack	75004117	6	9	12

Mattorna säljs i förpackningar om 2. Antal mattor för en plattform avser inte denna förpackning, utan en enstaka matta.

Tabell 46: Tillgängliga klistermattor för Solaris 4000-plattform

Tillbehör

### För Solaris 2000 I / 2000 R-plattformar

	Artikel- nummer	Universal
200 x 200 mm matta med hög vidhäftning, 2-pack	75004126	1
280 x 170 mm matta med hög vidhäftning, 2-pack	75004127	2
200 x 200 mm matta med låg vidhäftning, 2-pack	75004111	1
280 x 170 mm matta med låg vidhäftning, 2-pack	75004117	2

Mattorna säljs i förpackningar om 2. Antal mattor för en plattform avser inte denna förpackning, utan en enstaka matta.

Tabell 47: Tillgängliga klistermattor för Solaris 2000 I/2000 R-plattform

#### För Solaris 4000 I / 4000 R-plattformar

	Artikel- nummer	Universal
		18 x 18
200 x 200 mm matta med hög vidhäftning, 2-pack	75004126	4
280 x 170 mm matta med hög vidhäftning, 2-pack	75004127	2
200 x 200 mm matta med låg vidhäftning, 2-pack	75004111	4
280 x 170 mm matta med låg vidhäftning, 2-pack	75004117	2

Mattorna säljs i förpackningar om 2. Antal mattor för en plattform avser inte denna förpackning, utan en enstaka matta.

Tabell 48: Tillgängliga klistermattor för Solaris 4000 I/4000 R-plattform

Tillbehör

### 1. 2. 12. Allmänna tillbehör

I. 2. 12. Allmänna tillbehör								
	Artikelnummer							
Roller för klistermatta	75004124							
Skruvmejsel för tillbehör	75004131							
Gasfördelare för Solaris 2000 I/2000 R-plattform	SK2000-8GM							
Gasfördelare för Solaris 4000 I/4000 R-plattform	SK4000-8GM							
Klämma, Reservdelssats (Skruvar)	SK0010							
Reservdelssats för plattform för Solaris 2000, 2000 I/R och 4000 I/R (plattformsskruvar, verktyg)	SK0100							
Plattform reservdelssats för SK4000 (plattformsskruvar, verktyg)	SK0101							

Tabell 49: Tillgängliga allmänna tillbehör

### 1.3. Direktiv och standarder

Region	Direktiv	Standarder			
Europa	2006/42/EG	EN 61010-1			
	Maskindirektivet	IEC 61010-2-051			
	2011/65/EU RoHS	EN 61326-1 klass B			
	Direktiv för användning av farliga	EN ISO 14971			
	ämnen i elektronisk och elektrisk utrustning	EN ISO 9001			
	Skyddsmål:	Endast Solaris 2000 I/2000 R/			
	2014/35/EU Lågspänning	<u>4000 I/4000 R:</u> IEC 61010 2 010			
	2014/30/EG Elektromagnetisk kompatibilitet (EMC)				
Nordamerika		ANSI/UL 61010-1			
		IEC 61010-2-051			
		IEC 61326-1 klass B			
		CFR 47 FCC 15 EMC			
		EN ISO 14971			
		EN ISO 9001			
		Endast Solaris 2000 I/2000 R/			
		<u>4000 I/4000 R:</u> IEC 61010-2-010			
Japan		IEC 61010-1			
		IEC 61010-2-051			
		IEC 61326-1 klass B			
		EN ISO 14971			
		EN ISO 9001			
		Endast Solaris 2000 I/2000 R/ 4000 I/4000 R: IEC 61010-2-010			

Tabell 50: Direktiv och standarder

OBS! Denna utrustning har testats och befunnits överensstämma med gränserna för en digital enhet av klass A, i enlighet med del 15 av FCC-reglerna. Dessa gränser är utformade för att ge rimligt skydd mot skadliga störningar när utrustningen används i en kommersiell miljö. Denna utrustning genererar, använder och kan utstråla radiofrekvent energi och kan, om den inte installeras och används i enlighet med bruksanvisningen, orsaka skadliga störningar på radiokommunikationer. Användning av utrustningen i ett bostadsområde medför sannolikt störningar, varvid användaren måste korrigera störningen på egen bekostnad.

### 1.4. Produktöversikt

### 1.4.1. Solaris 2000



① Plattform, monteringsplatta; ② Grafiskt användargränssnitt (GUI); ③ Nätströmbrytare;
 ④ Nätanslutning; ⑤ Säkring; ⑥ Ethernet-port; ⑦ USB-port;

Bild 1: Översikt Solaris 2000

#### 1.4.2. Solaris 4000



① Plattform, monteringsplatta; ② Grafiskt användargränssnitt (GUI); ③ Nätströmbrytare;

④ Nätanslutning; ⑤ Säkring (återställningsbar); ⑥ Ethernet-port; ⑦ USB-port;

Bild 2: Översikt Solaris 4000

Produktöversikt

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



① Plattform, monteringsplatta; ② Grafiskt användargränssnitt (GUI); ③ Nätströmbrytare;

④ Tillbehörsport; ⑤ Nätanslutning; ⑥ Säkring (återställningsbar); ⑦ Ethernet-port;

⑧ USB-port; ⑨ Lock

Bild 3: Översikt Solaris 2000 I / 2000 R

### 1.4.4. Solaris 4000 I / 4000 R



① Plattform, monteringsplatta; ② Grafiskt användargränssnitt (GUI); ③ Nätströmbrytare;
 ④ Tillbehörsport; ⑤ Nätanslutning; ⑥ Säkring (återställningsbar); ⑦ Ethernet-port;
 ⑧ USB-port; ⑨ Lock

Bild 4: Översikt Solaris 4000 I / 4000 R

### 1.4.5. Anslutningar

### Öppna skakapparater



① Nätanslutning; ② Säkring; ③ Ethernet-port

Bild 5: Bakifrån (vänster: Solaris 4000, höger: Solaris 2000)

# Inkuberade skakapparater (uppvärmda) och kylda skakapparater (kylda och uppvärmda)



① Nätanslutning; ② Säkring; ③ Ethernet-port; ④ Dräneringsrör

Bild 6: Bakifrån (vänster: Solaris 2000 I / 2000 R, höger: Solaris 4000 I / 4000 R)

### <u>Nätanslutning</u>

Skakapparaten kräver strömförsörjning enligt tabellen nedan. Strömförsörjningskablar medföljer.

Artikel- nummer	Skakapparat	Specifikationer
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

Bild 7: Strömförsörjning för skakapparater

VARNING Felaktig nätanslutning eller nätkontakt medför skador på utrustningen. Se till att skakapparaten är ansluten endast till nätuttag som är korrekt jordat enligt gällande elsäkerhetsbestämmelser. Använd inte skakapparaten med en skadad eller otillräckligt klassad strömförsörjningskabel.

Stickproppen måste hela tiden vara fritt tillgänglig.

För att eliminera risken för elektriska stötar, se till att ytan runt skakapparaten är torr. I händelse av oavsiktligt spill eller vätskestänk, koppla bort skakapparaten från spänningskällan, torka upp spillet och eliminera hälsorelaterade eller biologiska risker innan du fortsätter.

Koppla bort skakapparaten från strömkällan när den inte används.

#### <u>Säkring</u>

Säkringen kan återställas om den har utlösts till följd av överström. När säkringen har återställts till sitt ursprungliga läge fungerar skakapparaten igen.

#### Ethernet

Skakapparaten har en RJ45 Ethernet-port som kan användas för anslutning till ett lokalt nätverk. Anslut endast utrustning som efterlever standarden IEC 60950-1 till Ethernet-porten. Ethernet-porten kan användas för framtida uppdatering av programvaran.

#### <u>Tillbehörsportar</u>

Temperaturstyrda modeller av skakapparaten är utrustade med två tillbehörsportar. Dessa är normalt förslutna med avtagbara lock. Locken kan tas bort om skakapparaten ska kompletteras med ytterligare instrument, exempelvis en extra temperaturgivare eller kabel- eller rörgenomföringar såsom gasfördelare.

### <u>USB</u>

Skakapparaten har två USB-A 2.0-portar för anslutning av USB-minne. Använd endast utrustning som överensstämmer med IEC 60950-1-standarden med USB-portarna.

#### Inbyggt överhettningsskydd

Temperaturstyrda modeller av skakapparaten är utrustade med ett manuellt återställningsbart termostat intill värmeelementen. Termostatet stänger av värmeelementen för att förhindra alltför höga temperaturer i skakapparatens kammare i händelse av att luftcirkulationsfläktarna skulle sluta fungera.

Termostatet måste återställas manuellt innan skakapparaten kan användas igen. Detta kan göras endast av Thermo Fisher Scientifics kundtjänst.

## 2. Transport och uppställning



Det är ditt ansvar att se till att skakapparaten ställs upp korrekt.

#### OBS

Transportförpackningen ska inspekteras vid leverans. Undersök leveransen noga och kontrollera eventuella fraktskador före uppackning. Om någon skada upptäcks ska transportföretaget specificera och underteckna skadan på din kopia av kvittot.

Öppna emballaget och kontrollera att alla delar (se "Tabell 51: Leveransomfattning") är inkluderade innan förpackningen kasseras. Om någon skada upptäcks efter uppackning: Rapportera det till transportören och begär en skadeinspektion.

Viktigt: Underlåtenhet att begära en inspektion av skada inom några dagar efter mottagandet av transporten befriar transportören från allt ansvar för skador. Det är ditt ansvar att kräva en skadeinspektion.

### 2.1. Uppackning

Använd packlistan vid uppackning för att verifiera att leveransen är komplett. Kassera inte förpackningsmaterial förrän du fastställt att alla delar finns med.

### Leveransomfattning

Artikel	Antal
Skakapparat	1
Nätkabel	1
Universalplattform	1*
Skruvar till plattform	
– Solaris 4000	4*
– Övriga modeller	3*
Tryckt engelskspråkig	1
bruksanvisning	
Bruksanvisning på USB	1
Verktyg (skruvnyckel med T-handtag)	1
* 2000 I-, 2000 R-, 4000 I- och 4000 R-mo	dellerna levereras med
förinstallerad plattform. Varken plattform elle	er tillhörande skruvar
medfölier som separata delar	

Tabell 51: Leveransomfattning

Om några objekt saknas, kontakta Thermo Fisher Scientific.

### 2.2. Uppställningsplats

SE UPP Skyddet kan försämras på grund av minskad stabilitet när plasten utsätts för ultraviolett strålning. Utsätt inte skakapparat eller plasttillbehör för direkt solljus eller andra källor av ultraviolett strålning.

Ställ skakapparaten på ett horisontellt avvägt bord eller bänk som kan bära skakapparatens vikt med tillbehör och prover under drift. Placera skakapparaten i närheten av ett eluttag som motsvarar kraven på märkskylten. Ge utrymme runt enheten för fri luftkonvektion och tillbehör. Gör arbetsområdet bekvämt för användaren.

Beakta följande krav för att sätta upp enheten:

- Minst 8 cm (3 in) fritt utrymme på plattformens alla fyra sidor krävs för att säkerställa tillräckligt med plats för rörliga delar, och för att undvika klämskador eller skador på närliggande utrustning. Den öppna skakapparatens plattform kan skjuta ut utanför höljet.
- Skakapparater ger upphov till vibrationer. Förvara inte känslig utrustning eller farliga föremål eller substanser inom säkerhetsområdet.



Bild 8: Öppen skakapparat måste ha minst 8 cm (3 in) fritt utrymme på alla fyra sidor

Följande utrymmeskrav måste vara uppfyllda för temperaturstyrda skakapparater för att säkerställa betryggande ventilation och optimala prestanda:

- För de inkuberade (uppvärmda) skakapparaterna Solaris 2000 I och 4000 I krävs minst 8 cm (3 in) fritt utrymme på alla fyra sidorna, (till vänster i Bild 9).
- För de kylda skakapparaterna (kylda och uppvärmda) Solaris 2000 R och 4000 R är 8 cm (3 in) fritt utrymme på alla fyra sidorna tillräckligt om utrymmet ovanför locket är fritt (till vänster i Bild 9).

 Om utrymmet ovanför de kylda skakapparaterna (kylda och uppvärmda) Solaris 2000 R och 4000 R är blockerat måste utrymmet runt höljet ökas till 30 cm (12 in) på alla fyra sidorna (till höger i Bild 9).



**Bild 9:** Installationskrav för inkuberade (uppvärmda) skakapparater (vänster) och kylda skakapparater (kylda och uppvärmda) (höger); kylda skakapparater (kylda och uppvärmda) med fritt utrymme ovanför illustreras till vänster i bilden.

VARNING Farliga ämnen måste hållas utanför denna zon under skakningen.

Ytan där apparaten ställts upp måste:

- » vara ren,
- » vara stabil, solid, stel och fri från resonans,
- » vara lämplig för horisontell uppställning av skakapparaten,
- » hålla skakmaskinens vikt.
- Skakapparaten får endast användas inomhus.
- Skakapparaten får inte utsättas för värme och starkt solljus.
- Dippställningsplatsen måste vara väl ventilerad hela tiden.
- Stickproppen måste hela tiden vara fritt tillgänglig.

### 2.3. Transportera

SE UPP Att tappa skakapparaten kan vålla fysiska skador. Lyft alltid skakapparaten från båda sidorna. Lyft aldrig skakapparaten i frontpanelen eller i en monterad plattform. Lyft aldrig en temperaturstyrd skakapparat i det transparenta lockets handtag.

SE UPP Säker användning av skakapparaten försämras om vatten har kondenserats på höljets insida. Detta kan hända om skakapparaten flyttas från en kallare till en varmare omgivning med fuktigare luft, till exempel från ett lagerutrymme eller motsvarande till laboratoriet. Låt utrustningen torka ut i minst 2 timmar före drift.

OBS Med öppna skakapparater ska plattform, last och alla tillbehör avlägsnas innan skakapparaten flyttas. Om du inte tar bort plattformen kan du skada plattformens monteringsplatta eller skakmekanismen. Med temperaturstyrda skakapparater är det tillräckligt att avlägsna lasten från plattformen.

**OBS** Stötar och slag kan skada enheten. Transportera skakapparaten i upprätt läge och om möjligt i originalförpackningen.

### Hantering

Tänk på följande vid hantering av skakapparaten:

- Lyft skakapparaten i dess sidor; aldrig i dess fram- och baksida.
- Minst två personer ska hjälpas åt att bära skakapparaten. Beroende på vikt (se "Specifikationer" på sidan 11) kan fler än två personer krävas.





Bild 10: Hur man bär en skakapparat

VARNING Lyft alltid skakapparaten från båda sidorna. Lyft aldrig skakapparaten i dess fram- och eller baksida. Skakapparaten är tung (se "Specifikationer" på sidan 11). Minst två personer krävs för att lyfta och bära skakapparaten. SE UPP Lyft eller bär aldrig en temperaturstyrd skakapparat i det transparenta lockets handtag. Handtaget riskerar att brytas sönder så att skakapparaten faller ned och skadar egendom eller människor.

### 2.4. Avvägning

Skakapparaten kan inte avvägas av sig själv. Skakapparaten kräver en avvägd horisontell yta för korrekt inställning.

OBS Placera inte mellanlägg av något slag mellan skakapparatens fötter och underlaget för att avväga den.

### 2.5. Nätanslutning

Skakapparaten kräver strömförsörjning enligt tabellen nedan. Strömförsörjningskablar medföljer. Se Bild 7 på sidan 47 för detaljer.

VARNING Felaktig nätanslutning eller nätkontakt medför skador på utrustningen. Se till att skakapparaten är inkopplad endast i uttag som är ordentligt jordade. Använd inte skakapparaten med en skadad eller otillräckligt klassad strömförsörjningskabel.

**OBS** Elektromagnetisk strålning kan ge upphov till störningar på displayen. Detta medför ingen risk för skada, begränsad funktion eller annan oönskad inskränkning. Undvik störningar från elektromagnetisk strålning genom att inte placera mobiltelefoner eller annan elektronisk utrustning i skakapparatens omedelbara närhet. Använd inte skakapparaten samtidigt med andra effektkrävande apparater som är anslutna till samma säkringsgrupp. Anslut inte flera apparater till ett och samma grenuttag.

Gör så här för att ansluta skakapparaten till strömförsörjningen:

- 1. Stäng av nätströmbrytaren som finns på höger sida.
- 2. Kontrollera att kablaget uppfyller gällande säkerhetsbestämmelser.
- 3. Försäkra dig om att centrifugens nätspänning och frekvens enligt märkskylten överensstämmer med den aktuella nätanslutningen.

Stickproppen måste hela tiden vara fritt tillgänglig.

För att eliminera risken för elektriska stötar, se till att ytan runt skakapparaten är torr. I händelse av oavsiktligt spill eller vätskestänk, koppla bort skakapparaten från spänningskällan och torka upp spillet innan du fortsätter.

Koppla bort skakapparaten från strömkällan när den inte används.

### 2.6. Första start

Utför följande steg innan du använder enheten:

1. Slå på strömmen till enheten. Thermo scientific-logotypen visas. Tryck **Start Setup**.



Bild 11: Skärmen vid första start

2. Välj önskat språk på skärmen Språk. Tryck Nästa.



Bild 12: Första start - språk

 Om så önskas kan ett namn på enheten anges i dialogrutan Enhetsnamn. Tryck Nästa.

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

Bild 13: Första start - enhetsnamn

 Om så önskas kan ort och land anges i textfältet. När du har skrivit de tre första tecknen visas en lista över matchande förslag i textfältet Ort, Land. Välj ett passande förslag och tryck på Nästa.

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

Bild 14: Första start - region

5. Välj önskat datumformat. Tryck Nästa.

Set Date Dis	play Format
	I/DD/YYYY
O DD/M	IMM/YYYY
O YYYY	/MMM/DD
Back	Next

Bild 15: Första start - ange visningsformat för datum

6. Tryck på respektive upp- eller nedpil eller dra med fingret för att ställa in aktuellt datum. Tryck Nästa.



Bild 16: Första start – ställ in datum

 Välj önskat tidsformat och tryck på respektive upp- eller nedpil eller dra med fingret för att ställa in aktuell tid. Tryck Nästa.



Bild 17: Första start – ställ in tid

 Endast inkuberade skakapparater (uppvärmda) och kylda skakapparater (kylda och uppvärmda): Välj önskad temperaturenhet. Tryck Nästa.

Temperature Units								
⊙* <b>●</b>								
○ 'F								
Back Next								
Dack								

Bild 18: Första start – välj temperaturenhet

 Endast inkuberade skakapparater (uppvärmda) och kylda skakapparater (kylda och uppvärmda): Tryck på den blå pilen för att ställa in larm om hög temperatur och tröskelvärden för nödavstängning. Tryck Nästa.



Bild 19: Första start – ställ in tröskelvärden för larm om hög respektive låg temperatur

10. Upprepa processen för tröskelvärden för larm om låg temperatur. Tryck på Nästa för att fortsätta.

11. Om du vill fortsätta utan att definiera ett administratörslösenord väljer du **Öppet läge** och trycker på **Nästa** för att fortsätta.

Om du vill definiera ett administratörslösenord trycker du på Skyddat läge.



Bild 20: Första start - välj behörighetsläge

12. En lösenordsskärm visas; ange önskat administratörslösenord via tangentbordet på skärmen och tryck på **Nästa**.



Bild 21: Första start – välj behörighetsläge

- 13. På nästa lösenordsskärm som visas anger du administratörslösenordet.
- 14. En tredje lösenordsskärm visas, där du uppmanas att bekräfta det nya administratörslösenordet genom att ange det ännu en gång.

	Admin Passcode					Admin Passcode									
If you	do not wis fields	h to change below blanl	the pa c and ta	sscode ap the N	at this ti ext butto	ime, leav on.	e the		Ρ	lease re-ent	er the ne	ew pass	code.		
	*	*	*		*	*			*	*	*	-	e.	*	
		7	8	9						7	8	9			
		4	5	6						4	5	6			
		1	2	3						1	2	3			
		0	CI	ear						0	Cle	ar			

Bild 22: Första start – ange det nya administratörslösenordet och bekräfta

15. Tryck på Nästa på bekräftelseskärmen för att fortsätta.

lode					
fully. Please keep the new fe location as you will need it ettings.					
If the passcode is forgotten you will need to contact Customer Service to reset.					
Next					
	ode fully. Please keep the new e location as you will need it titings. ill need to contact Customer reset. Next				

Bild 23: Första start – bekräfta ändring av administratörslösenordet

16. En skärm med allmänna instruktioner för fysisk installation av skakapparaten visas. Tryck Nästa.

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			

Bild 24: Första start - Installationsanvisningar

17. Skärmen för slutförd installation visas. Tryck på Nästa för att slutföra.

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

Bild 25: Första start – utför kalibrering eller avsluta konfigureringen

**OBS** Med inkuberade skakapparater (uppvärmda) och kylda skakapparater (kylda och uppvärmda) kan du trycka på **Kalibrera** och utföra proceduren som beskrivs i kapitel "4. 6. Temperaturkalibrering" på sidan 126.

### 2.7. Förvaring



SF UPP

När du tar bort skakapparaten och tillbehören från användning, rengör och vid behov desinficera eller dekontaminera hela systemet. Lämna inte skakapparaten och tillbehör i ett odefinierat tillstånd av förorening. Om du är osäker på processen, kontakta Thermo Fisher Scientific kundtjänst ("Rengöring" på sidan 123, "Desinficering" på sidan 124 och "Dekontaminering" på sidan 124).

- Innan skakapparaten och tillbehören ställs undan skall de rengöras och vid behov desinficeras och dekontamineras.
- Skakapparaten och tillbehören måste vara helt torra före förvaring.
- Förvara skakapparaten på en ren, dammfri plats.
- Se till att skakapparaten står på fötterna.
- Förvara inte skakapparaten i direkt solljus.

### 2.8. Leverans



SE UPP

Innan du transporterar skakapparaten och tillbehören måste du rengöra och vid behov desinficera eller dekontaminera hela systemet. Lämna inte skakapparaten och tillbehör i ett odefinierat tillstånd av förorening. Om du är osäker på processen, kontakta Thermo Fisher Scientific kundtjänst ("Rengöring" på sidan 123, "Desinficering" på sidan 124 och "Dekontaminering" på sidan 124).

Innan du transporterar skakapparaten:

- Skakapparaten måste vara ren och dekontaminerad.
- Dekontaminationen måste verifieras med ett dekontamineringscertifikat. Ett dekontaminerinsintyg kan fås från Thermo Fisher Scientific kundtjänst.

# 3. Drift

### 3.1. Ström på/av

Tryck på strömbrytaren på höger sida för att starta (I) eller stänga av (0) skakapparaten.

Under startsekvensen visas Thermo Scientific-logotypen på pekskärmen.

När den är klar, visar pekskärmen aktuell status för skakapparaten.

### 3.2. Grafiskt användargränssnitt

Startskärmen är standardskärmen i skakapparatens grafiska användargränssnitt. Den utgör startpunkten för användning av skakapparaten.

På startskärmen kan du:

- Ställa in grundläggande driftinställningar såsom skakningshastighet, tid och temperatur (endast temperaturstyrda skakapparater).
- Starta och stoppa skakapparaten.
- Visa statusinformation och hantera larm och varningar.
- Gå till andra skärmar med statusinformation och inställningar.

Skärmens utseende skiljer sig något mellan öppna och temperaturstyrda skakmodeller.



Ett exempel på startskärm för öppen skakapparat visas i Bild 26.

Bild 26: Startskärm för öppen skakapparat

Startskärmen för en temperaturstyrd skakapparat har extra funktioner för temperaturstyrning (Bild 27).



Bild 27: Startskärm för temperaturstyrd skakapparat

### 3. 2. 1. Grundläggande driftinställningar

Från startskärmen når du med en knapptryckning inställningsskärmarna för skakapparatens huvudsakliga driftinställningar.

#### Hastighetsinställning

1. Tryck på fältet **Faktisk hastighet** (**(i** Bild 26 eller **()** i Bild 27) för att öppna skärmen **Hastighetsinställningar** (Bild 28).



Bild 28: Hastighetsinställningar

- 2. Tryck på respektive uppil eller nedpil för att ställa in önskad hastighet.Du kan också ändra hastigheten genom att dra värdena uppåt eller nedåt med fingret.
- 3. Tryck Spara.

**OBS** Om ett meddelande visas att hastighetsinställningen är ogiltig är inställningen utanför skakapparatens möjliga drifthastighetsområde (kapitlet "Tekniska data" på sidan 11). Korrigera inställningen och fortsätt.

4. Bekräfta dialogrutan Ändringarna har sparats för att återgå till huvudskärmen.

### Beräkna omkretsbana: Beräkna approximativ hastighet från inställningarna för äldre utrustning

I dialogrutan Hastighetsinställningar kan du trycka på **Beräkna hastighet** för att visa skärmen **Beräkna omkretsbana.** Denna funktion är praktisk vid övergång från en skakapparat med en annan mekanisk konstruktion till en av Solaris-skakapparaterna som beskrivs i den här bruksanvisningen. Denna kalkylator använder omkretsbanans differens för att beräkna en ungefärlig hastighet som krävs för att uppnå ett motsvarande resultat.

OBS Använd inte det resultat som erhålls med funktionen Beräkna omkretsbana för dina prover utan att först bekräfta resultatet via en serie tester. Celler kan växa snabbare eller långsammare, utsöndra annorlunda proteiner eller skadas eller dö till följd av de ökade fysiska påfrestningar som kan uppkomma när en annorlunda omkretsbana används.

1. Välj måttenhet i millimeter eller tum.



Bild 29: Beräkna omkretsbana

 Välj storlek på föregående omkretsbana. Tryck **Beräkna**. Följande skärm för hastighetsberäkning visas.



Bild 30: Beräknad omkretsbana

#### 3. Tryck OK.

- Du kan beräkna den föreslagna hastigheten för en anpassad storlek på omkretsbanan. Om du vill göra en beräkning för en anpassad omkretsbana väljer du Annan.
- 5. Ange omkretsbanans storlek.



Bild 31: Beräkna omkretsbanan för anpassad storlek

6. Tryck Beräkna.

Fönstret Beräkna hastighet visas med hastigheten för den angivna storleken.

#### Ställa in körtid

Skakapparaten kan köras i läget Kontinuerligt eller Tidsinställt. I läget Kontinuerligt måste skakapparaten stoppas manuellt. I läget Tidsinställt stoppar skakapparaten när den inställda tiden har löpt ut. Timern kan visas på olika sätt:

- Förfluten tid: den tid som skakapparaten har körts sedan du tryckte på Start.
- Återstående tid: den tid som skakapparaten kommer att köra tills den inställda tiden har löpt ut.
- 1. Tryck på fältet **Tid** (⑤ i Bild 26 och Bild 27) för att öppna skärmen Tidsinställt (Bild 32).
- 2. Välj Tidsinställt.



Bild 32: Välj Tidsinställt

- 3. Ange timmar och minuter i rutorna Ställ in tid.
- 4. Tryck Spara.
- 5. Bekräfta dialogrutan Ändringarna har sparats för att återgå till huvudskärmen.

#### Ställa in temperatur

Med temperaturstyrda modeller kan du när som helst ställa in en temperatur för den aktuella uppgiften i fältet **Temperaturbörvärde**.

- 1. Tryck på fältet **Temperaturbörvärde** (6 Bild 27) för att visa skärmen **Temperaturbörvärde** (Bild 33).
- Tryck på respektive upp- eller nedpil för att förinställa önskad standardtemperatur för skakapparaten.Du kan ändra temperaturen även genom att dra värdena uppåt eller nedåt med fingret.



Bild 33: Ställ in temperaturbörvärde: Skärm med kalibreringsvärde

3. Tryck Spara.

**OBS** Om ett meddelande visas att temperaturbörvärdet är ogiltigt är inställningen utanför skakapparatens möjliga drifttemperaturområde (kapitlet "Tekniska data" på sidan 11. Korrigera inställningen och fortsätt.

**OBS** Om ett meddelande visas att den valda temperaturen är ogiltig på grund av den aktuella omgivningstemperaturen är omgivningstemperaturen utanför skakapparatens möjliga drifttemperaturområde (kapitlet "Tekniska data" på sidan 11). Tryck på **OK** för att bekräfta varningen och fortsätt, eller ställ in en annan temperatur.

OBS På skärmen **Temperaturbörvärde** visas den kalibrerade temperaturen och avvikelsen som justerades vid kalibreringen (nederst till vänster i Bild 33). Kalibreringsprocessen beskrivs i "Temperaturkalibrering" på sidan 126.

4. Bekräfta dialogrutan Ändringarna har sparats för att återgå till huvudskärmen.

OBS Temperaturstyrda skakapparater beter sig på ett specifikt sätt när de uppnår inställd temperatur, så kallat temperatursvängning. Temperatursvängning innebär att den inställda kammartemperaturen först överskrids till en något högre (eller lägre) temperatur för att därefter justeras till inställd temperatur. När den inställda temperaturen uppnås visas inte denna temperatursvängning på pekskärmen. Istället visas den stigande (eller fallande) kammartemperaturen tills inställd temperatur uppnås.

#### Starta och stoppa skakapparaten

1. Starta skakapparaten genom att trycka på Start.

Startknappen förvandlas till Stoppknapp.



Bild 34: Start- och stoppknapp

2. Stoppa skakapparaten genom att trycka på **Stopp**.

### 3. 2. 2. Status

När skakapparaten fungerar normalt visas en grön hjärtikon i området Hälsostatus



och statistik (⑦ i Bild 26 och ⑧ i Bild 27) på skärmen). Om du trycker på den gröna hjärtikonen öppnas statusskärmen. Statusskärmen innehåller driftinformation om skakapparaten, inklusive antal skaktimmar och

drifttimmar totalt och för den senaste sessionen. För temperaturstyrda skakapparater visas dessutom antal kylnings- eller uppvärmningstimmar.



**Bild 35:** Statistisk information: Öppen skakapparat (vänster) och temperaturstyrd skakapparat (höger)

### Meddelande

När ett meddelande visas, visas ett gult fält överst på skärmen. Dessutom avges en ljudsignal. Det gula meddelandefältet försvinner när du rullar innehållet i

meddelandelistan två gånger. Endast den gula triangeln indikerar att det finns ett eller flera meddelanden. I den triangelformade ikonen finns en blå cirkel med en vit ram som indikerar antalet aktiva meddelanden. Genom att



cirkel med en vit ram som indikerar antalet aktiva meddelanden. Genom att trycka på varningstriangeln i området Hälsostatus och statistik (⑦ i Bild 26 och ⑧ i Bild 27) visas en skärm med alla aktiva meddelanden. Det senaste meddelandet visas utökat så att du kan se alla detaljer (Bild 36). Du kan bläddra i listan och tryck på ett listobjekt för att expandera det och läsa mer.



Bild 36: Lista med meddelanden

Välj det aktiva meddelandet genom att trycka på kryssrutan intill meddelandet. Tryck på **Kvittera** för att ta bort meddelandet från listan. När alla meddelanden har tagits bort återgår statusikonen till den gröna hjärtikonen.

#### Larm

När ett larm utfärdas stannar skakapparaten omedelbart för att undvika skador på



proverna och/eller själva enheten. Du måste kvittera larmet på pekskärmen innan driften kan fortsätta.

När ett larm utfärdas visar pekskärmen ett rött fält ovanpå den aktuella skärmen. I området Hälsostatus och statistik (⑦ i Bild 26 och ⑧ i Bild 27) visas en röd larmklocka. Dessutom avges en konstant ljudsignal.

Nedanför det röda larmfältet visas en sammanfattning av det aktuella larmet. Med snooze-knappen till höger kan du tillfälligt stänga av alarmet. Om larmtillståndet inte blir löst inom snooze-perioden, startar det akustiska larmet igen. Varaktigheten för snooze-perioden kan väljas i inställningarna. Detta förklaras i avsnittet "Larm och meddelanden" på sidan 69.

Genom att trycka på larmklockan i området Hälsostatus och statistik (⑦ i Bild 26 och ⑧ i Bild 27) visas en skärm med fullständig information om det aktiva larmet (Bild 37). Du kan bläddra i listan och tryck på ett listobjekt för att expandera det och läsa mer.



Bild 37: Larmlista

Välj det aktiva larmet genom att trycka på kryssrutan intill larmet. Tryck på **Kvittera** för att ta bort larmet från listan när larmorsaken har åtgärdats. När alla larm har tagits bort återgår statusikonen till den gröna hjärtikonen.

### Fel

När ett fel inträffar visas ett felmeddelande och skakapparaten stoppar omedelbart för att förhindra skador på proverna och/eller enheten. Skärmen blir röd och skakapparaten kan inte längre aktiveras. Ett felmeddelande och en felkod visas; se exempel i Bild 38.



Bild 38: Skärm med felmeddelande

Gör så här för att återuppta driften:

- 1. Notera felkoden som visas på skärmen.
- 2. Tryck på Kvittera för att stänga av larmsignalen.
- 3. Starta om skakapparaten genom att stänga av den och slå på den igen.
- 4. Om felmeddelandet fortfarande visas, kontakta Thermo Fisher Scientifics kundtjänst och uppge felkoden som visas på skärmen.

### 3. 2. 3. Inställningar

Det andra alternativet i navigeringsfältet är ikonen Inställningar. När du trycker på **Inställningar** visas nedanstående skärm. Inställningsskärmen innehåller ytterligare knappar, men alla knappar kan inte visas samtidigt på skärmen. Tryck på upp- eller nedpilen för att visa återstående knappar (till höger i Bild 39).



Bild 39: Skärmen Inställningar

### Larm och meddelanden

På skärmen Larm och meddelanden kan du välja hur och när larm och meddelanden ska utlösas.



Bild 40: Inställningar för larm och meddelanden

**Volym**: Med den här funktionen kan du ändra meddelandesignalens ljudvolym för ett larm eller meddelande som utlöses. Dra skjutreglaget åt vänster för att sänka, eller åt höger för att höja, volymen och tryck på **Spara**. Denna volyminställning påverkar både larm och meddelanden.



Bild 41: Ställa in ljudvolymen för larm och meddelanden

Larm hög temperatur/larm låg temperatur (endast temperaturstyrda skakapparater): Med dessa två alternativ kan du ställa in ett larmtröskelvärde för hög och låg temperatur i förhållande till temperaturbörvärdet ("Ställa in temperatur" på sidan 64). När temperaturen i skakapparatens kammare underskrider tröskelvärdet för låg temperatur, eller överskrider tröskelvärdet för hög temperatur, utlöses ett larm. Välj ett temperaturtröskelvärde och tryck på **Spara**.



Bild 42: Ställa in larm för hög och låg temperatur

OBS Om du ändrar temperaturbörvärdet justeras larmtröskelvärdena i motsvarande grad.

Avstängning vid hög temperatur/avstängning vid låg temperatur (endast temperaturstyrda skakapparater): Med dessa två alternativ kan du ställa in ett avstängningströskelvärde för hög och låg temperatur i förhållande till temperaturbörvärdet ("Ställa in temperatur" på sidan 64). När temperaturen i skakapparatens kammare underskrider tröskelvärdet för låg temperatur, eller överskrider tröskelvärdet för hög temperatur, stängs skakapparaten av automatiskt för att skydda proverna i skakapparatens kammare. Välj ett tröskelvärde och tryck på Spara.

High Temperature S	+15 🔨	Low Temperature Sh	-15 🔨
	+20		-20
	+25		-25

Bild 43: Ställa in avstängningströskelvärde för hög och låg temperatur

OBS Om du ändrar temperaturbörvärdet justeras tröskelvärdena i motsvarande grad.

**Snooze Timeout**: Tidsintervallet för Snooze-timeout avgör hur lång tid ett larm är tyst när du trycker på Snooze i det röda fältet överst på huvudskärmen ("Larm" på sidan 67). Du kan ställa in snooze-timeout på 5 minuter, 10 minuter eller 15 minuter. Standardalternativet är 10 minuter. Tryck på **Spara** för att spara ändringarna.



Bild 44: Ställa in snooze-timer för larm

Inaktivera kalibreringsmeddelanden (endast temperaturstyrda skakapparater): Med dessa alternativ kan du inaktivera meddelanden som regelbundet påminner om att kalibrera skakapparatens temperaturmätning ("Temperaturkalibrering" på sidan 126). Du kan välja att inaktivera kalibreringspåminnelser helt, eller endast när ett program körs.



Bild 45: Inaktivera kalibreringspåminnelser

### <u>Behörighetskontroll</u>

På skärmen Behörighetskontroll kan du definiera ett administratörslösenord för funktioner med krav på begränsad åtkomst.

Som standard är skakapparaten inställd på Öppet läge, dvs. inget lösenord krävs för att använda den. Om skakapparaten ställs in på Skyddat läge måste alla användare som vill ändra skakapparatens inställningar ange administratörslösenordet. I skyddat läge visas en inloggningsknapp i skärmens övre högra hörn.

Följande jämförelse mellan Öppet läge och Skyddat läge visar när en lösenordsskyddad inloggning krävs.

Åtgärd	Lösenord krävs i Öppet läge	Lösenord krävs i Skyddat läge
Ställa in skakningshastighet, tid och temperaturbörvärden	Nej	Nej
Starta och stoppa skakapparaten	Nej	Nej
Beräkna omkretsbanor	Nej	Nej
Kvittera larm och meddelanden	Nej	Nej
Visa skärmen Filer och information	Nej	Nej
Visa hälsostatus	Nej	Nej
Visa och exportera händelselogg och diagram	Nej	Nej
Snooza larm	Nej	Nej
Köra program	Nej	Nej
Skapa, redigera och ta bort program	Nej	Ja
Ändra visningsinställningar	Nej	Ja
Ändra kontrollinställningar	Nej	Ja
Ändra inställningar för larm och meddelanden	Nej	Ja
Ändra behörighetsinställningar	Nej	Ja
Importera eller exportera program	Nej	Ja
Återställa till fabriksinställningarna	Nej	Ja
Utföra kalibrering	Nej	Ja
Installera uppdatering av inbyggd programvara	Ja	Ja

Tabell 52: Krav på lösenordsskyddad inloggning i Öppet och Säkert läge

#### Grafiskt användargränssnitt

I Bild 46 visas skärmen Behörighetskontroll i öppet och skyddat läge.

< 余	Shake Lab 1 11:22:24AM Access Control		< *	Shake Lab 1 11:22:24AM Access Control	Logout Admin
¢.	<ul> <li>Open</li> <li>Secure Will require passcode</li> </ul>	access to	۵.	Open Secure Will require pas	scode access to
	change certain param Change Admin Passcode	eters		change certain j	parameters
<b>.</b>	Save	()		Save	

Bild 46: Behörighetskontroll: Öppet läge (vänster) och skyddat läge (höger)

När du trycker på **Skyddat** och därefter på **Spara** måste du ange administratörslösenordet för att bekräfta ändringen. Alla enheter levereras med samma administratörslösenord förinställt på fabrik. Lösenordet återfinns i bruksanvisningen.

Även användare som försöker ändra en lösenordsskyddad inställning uppmanas att ange administratörslösenordet.

OBS Det förinställda administratörslösenordet är 00000.

Med knappen Ändra admin-lösenordet kan du ändra lösenordet. När du trycker på knappen Ändra admin-lösenordet, uppmanas du att först ange det befintliga lösenordet och därefter det nya lösenordet. Ange det nya lösenordet en gång till för att bekräfta. Meddelandet Nytt lösenord har sparats visas som bekräftelse.

#### Kalibrering

Via kalibreringsskärmen (endast temperaturstyrda skakapparater) visas en serie skärmar där skakapparaten kan kalibreras för exakt temperaturstyrning. Kalibreringsprocessen beskrivs i "Temperaturkalibrering" på sidan 126.
## Kontroller

Skärmen Kontroller innehåller alternativ för standardinställningar för skakapparatens drift.

Hastighetsinställningar: Hastigheten kan ställas in mellan 15 och 525 v/min. Dra hastighetsvärdena uppåt eller nedåt med fingret för att förinställa hastigheten och tryck på **Spara**. En utförlig beskrivning av hur du använder skärmen Hastighetsinställningar finns i avsnittet "Hastighetsinställning" på sidan 61 kapitel 3. 2. 1.

**Tidsläge:** Skakapparaten kan köras i läget Kontinuerligt eller Tidsinställt. I läget Kontinuerligt måste skakapparaten stoppas manuellt. I läget Tidsinställt stoppar skakapparaten när den inställda tiden har löpt ut. I läget Tidsinställt kan du ställa in en standardkörtid och välja om förfluten tid eller återstående tid ska visas när skakapparaten körs. En utförlig beskrivning av hur du använder skärmen Tidsinställt finns i avsnittet "Ställa in körtid" på sidan 63 kapitel 3. 2. 1.

**Temperaturbörvärde** (endast temperaturstyrda skakapparater): Med detta alternativ kan du ställa in en standardtemperatur som ska användas när skakapparaten startar. En utförlig beskrivning av hur du använder skärmen Temperaturbörvärde finns i avsnittet "Ställa in temperatur" på sidan 64 kapitel 3. 2. 1.

Automatisk omstart: Den här funktionen startar om enheten efter ett strömavbrott vid normal start/stopp, programkörning eller temperaturkalibrering. Om den automatiska omstarten är inställd på **Nej** kommer enheten inte att startas om efter ett strömavbrott.

### <u>Display</u>

Med displayinställningarna kan du ändra olika visningsalternativ.

Ljusstyrka: Om du vill justera ljusstyrkan för displayen använder du glidkontrollen eller +/--knapparna.



Bild 47: Ställa in skärmens ljusstyrka

**Språk**: Om du vill ändra visningsspråk trycker du på **Språk**. Välj önskat språk genom att dra med fingret över värdena och trycka på **Spara**.



Bild 48: Språkinställning

Måttenhet (endast temperaturstyrda skakapparater): Tryck på Måttenhet för att välja Celsius °C eller Fahrenheit °F för temperaturavläsningar på alla skärmar.



Bild 49: Ställa in måttenhet för temperatur

**Datum**: Ställ in datumet genom att trycka på knappen Datum. Dra värdena för år, månad och dag uppåt eller nedåt med fingret för att ställa in aktuellt datum och tryck på **Spara**.



Bild 50: Ställa in år, månad och dag

**Datumformat**: Ställ in datumformat genom att trycka på knappen **Datumformat**. Tryck på ett av alternativen för önskat datumformat (exempel: DD/MM/YYYY visas som 15 april 2015) och tryck på **Spara**.



Bild 51: Ställa in datumformat

Tid: Om du vill ställa in Tid och tidsformat trycker du på knappen Tid. Tryck på 12 tim eller 24 tim och dra värdena för timme, minut och AM/FM (endast 12-timmarsformat) med fingret och tryck på **Spara**.



Bild 52: Ställa in tid och 12-timmars- eller 24-timmarsformat

Viloläge: Tryck på Viloläge om skakapparaten ska växla till viloläge efter 15 minuters inaktivitet. I viloläge visas meddelandet **Tryck för att vakna** efter de 15 minuterna.



Bild 53: Aktivera viloläge

**Region**: Ange i vilket land skakapparaten används genom att trycka på **Land**. Tryck på textfältet Ort, Land och börja skriva de första tecknen i ortnamnet. När du har skrivit de tre första tecknen visas matchande förslag i textfältet. Välj en matchande ort eller fortsätt skriva hela namnet och tryck på **Spara** på tangentbordet på skärmen.



Bild 54: Välj ort och land

Enhetsnamn: Om du vill definiera eller ändra skakapparatens namn trycker du på Enhetsnamn. Tryck på textfältet Enhetsnamn och skriv ett namn. Tryck på Spara på tangentbordet på skärmen.

< ♠	Sh 11 Ur	Shake Lab 1 11:22:24AM Unit Name										< ♠	Shake Lab 1 11:22:24AM Unit Name	$\bigcirc$	
۵.	Cunit Name				Shake Lab 1							۵. «			
Q	W	E	R	Т	Y	U	Ţ	0	Ρ			<b></b>	Unit Name:	Shake Lab 1	
А	S	D	F	G	Н	J	К	L	Sa	ave					
1	Z	X	С	۷	В	N	М	@		1		+			
?!123				-			_	<b>.</b>				Save	()		

Bild 55: Namnge skakapparaten

Anpassningsmenyn: Tryck på Anpassningsmenyn för att anpassa de två nedersta ikonerna i navigeringsfältet. Dra önskad ikon från huvudskärmen till den ikon som du vill ersätta. Tryck på Spara för att bekräfta.

OBS Du kan när som helst trycka på Återställ menyn om du vill återställa navigeringsfältet till fabriksinställningarna.



Bild 56: Anpassa navigeringsfältet

### Filer och information

På den här skärmen visas serienummer, användargränssnitt med versionsnummer, styrenhetens inbyggda programvara med versionsnummer, parameterfil med versionsnummer samt aktuellt operativsystem.

Du kan trycka på **Fabriksåterställning** om du vill återställa skakapparaten till fabriksinställningarna. Du måste ange administratörslösenordet; alla gjorda inställningar raderas. Händelseloggen raderas däremot inte.



Bild 57: Filer och information

Rulla nedåt i listan för att visa knappen **Inbyggd programvara** (till höger i Bild 57). Du kan uppdatera skakapparatens inbyggda programvara genom att trycka på denna knapp. Utförliga instruktioner om installation av ny inbyggd programvara finns i "4. 7. Installera inbyggd programvara" på sidan 130.

OBS Ny inbyggd programvara kan erhållas endast av en auktoriserad servicetekniker.

### <u>Service</u>

Serviceinställningarna kan endast nås av auktoriserade servicetekniker.

# 3. 2. 4. Program

Tryck på ikonen **Program** i navigeringsfältet för att visa en programlista eller skapa ett nytt program. Du kan skapa, redigera, ta bort, importera och exportera program. I Bild 58 visas skärmen Program för en ny enhet (vänster) och en enhet som har använts under en tid (höger). En lista med program skapade av användare visas för den aktuella enheten.

<	Shake Lab 1 11:22:24AM Programs	Shake Lab 1 11:22:24AM Programs
		Soluble Solution
۳°	Tap the Create New button to create a program.	Fizzy Bubbles 💉 💿
		martini 🖍 💿
		🔔 🕨 Champagne 💉 👁 🗸
	🕂 Create New Adv. Settings 🚺 🚺 🚺	😑 🕂 Create New Adv. Settings 🚺 💦 🕕

Bild 58: Program

Vid varje post i listan visas en ögonikon 🧿 som visar skärmen Snabbvisa program med programdetaljer, och en pennikon 📝 som visar programmet för redigering.

#### Gör inställningar för alla program

Tryck på **Avanc. inställn.** på Program-skärmen för att visa skärmen **Avancerade inställningar**. På den här skärmen kan du styra vad som ska ske när ett program startas. Dessa inställningar används för alla program du skapar.



Bild 59: Göra programinställningar

- 1. Tryck på Avanc. inställn.
- 2. Välj ett alternativ för Visa timer för att välja hur programtimern ska visas:
  - » Förfluten: den tid som skakapparaten har körts sedan du tryckte på Start.
  - » Återstående: den tid som skakapparaten kommer att köra tills den inställda tiden har löpt ut.

 Om skakapparaten ska börja köra omedelbart när ett program startas väljer du Starta skakapparaten när program väljs. Som standard måste du trycka på Start för att starta ett program.

OBS Om **Starta skakapparaten när program väljs** ska fungera måste skakapparatens lock vara stängt. När locket stängs startar körningen automatiskt.

4. Tryck Spara.

#### Skapa program

Du kan skapa och lagra upp till 99 program.

- 1. Tryck på Skapa nytt.
- 2. Ange namnet på programmet.





3. Ange temperatur (endast temperaturstyrda skakapparater), hastighet samt körtid för programmet i timmar och minuter.

Om temperaturstyrning hos en temperaturstyrd skakapparat ska inaktiveras lämnar du fältet **Temp** tomt.

- Om du vill lägga till ett extra steg i programmet bläddrar du nedåt och trycker på Lägg till steg- knappen.
- 5. Tryck på **Spara** för att spara programmet.

#### Redigera program

- 1. Tryck på pennikonen 🖍 bredvid det program du vill redigera.
- 2. Redigera önskade fält. Tryck **Spara**. Programmet sparas med de nya ändringarna.
- Du kan lägga till ett steg genom att välja Lägg till steg längst ned på skärmen. Bläddra på skärmen för att se knappen Lägg till steg om det finns mer än 3 steg.

OBS Om ett larm eller felmeddelande utfärdas under en körning i flera steg, stoppar skakapparaten automatiskt. Om ett meddelande utfärdas, kommer skakningen att fortsätta.



Bild 61: Lägg till steg

### Radera program

- 1. Tryck på pennikonen bredvid det program du vill redigera.
- 2. Tryck Ta bort. Ett fönster visas där du ombeds bekräfta.





3. Tryck **Ta bort**. Det valda programmet raderas.

#### Köra ett program

- 1. Om du vill köra ett befintligt program väljer du det program du vill köra.
- 2. Om du vill se snabbvisning av programmet, tryck på ögonikonen () bredvid programmet.

3. Ett fönster för snabbvisning öppnas.



Bild 63: Snabbvisa program

4. Tryck på Kör för att ladda programmet på startskärmen.

Beroende på valda inställningar för **Automatisk körning** startar programmet omedelbart, eller när du trycker på **Start** på startskärmen.

5. När programmet är avslutat visas följande fönster. Tryck **OK**.



Bild 64: Programmet är avslutat

#### Importera program

Du kan importera program som skapats i en skakapparat till en annan. Ett USBminne måste anslutas till skakapparatens USB-port.

På följande skärm visas alla program på det aktuella USB-minnet.



Bild 65: Välj program för import

Markera de program du vill importera. Tryck Importera.

När programmen har importerats visas följande fönster. Tryck **OK**. Du kan nu ta bort USB-enheten.



Bild 66: Programimport klar

### Exportera program

Du kan exportera program från en skakapparat till en annan. Kontrollera att en USBenhet är ansluten.

Följande skärm visas om USB-enheten är ansluten:



Bild 67: Välj program för export

Markera de program du vill exportera. Tryck Exportera.

Följande fönster visas när exporten är slutförd. Tryck **OK**. Du kan nu ta bort USBenheten.



Bild 68: Programexport klar

## <u>Händelselogg</u>

Den tredje fliken på navigeringspanelen är händelseloggen som innehåller användaroch systemhändelser. Tryck på ikonen Händelselogg i navigeringsfältet för att visa skärmen Händelselogg (Bild 69).



Bild 69: Händelselogg

På den här skärmen visas en lista över de senaste händelserna, med datum- och tidsstämpel för varje händelse.

Du kan visa ytterligare information om en enskild händelse genom att välja händelsen i listan.

Händelser kan filtreras efter typ enligt följande kategorier:

- larm
- meddelanden
- ändrade inställningar
- start- och stopp-åtgärder
- programkörningar
- locköppningar (endast temperaturstyrda skakapparater)

När ett filter väljs ändras knappen Visa alla till höger till Filter PÅ (Bild 70).



Bild 70: Filtrerad händelselogg

#### Exportera händelseloggen

1. Välj de händelser som ska exporteras i listrutan **Exportera händelser**. Välj exportformat för loggen eller rapporten.



Bild 71: Välj händelselogg för export

2. Ett fördefinierat eller anpassat datumintervall kan väljas.

OBS Endast händelser från de sex senaste månaderna kan exporteras.

 Export kan göras med hjälp av en USB-enhet. Sätt i en USB-enhet för att lagra loggen eller rapporten. Tryck på knappen Exportera logg för att hämta loggen eller rapporten.



Bild 72: Sätt i USB-enheten för export

## <u>Diagram</u>

Hastighets- eller temperaturdata (endast temperaturstyrda skakapparater) kan visas över tid i diagramform. X-axeln visar tid och Y-axeln visar hastighet eller temperatur.

Med temperaturdiagram kan du plotta börvärde, faktisk temperatur och omgivningstemperatur över tid. Tryck på **Redigera** för att välja vilka av de tre kurvorna som ska visas i temperaturdiagrammet. Tidsintervallet kan omfatta en dag, sju dagar eller valfritt antal dagar för antingen hela dygnet eller valda timmar varje dag.

Tryck på **Uppdatera** för att uppdatera diaggramvisningen enligt valda inställningar och visa nya data som har registrerats sedan diagrammet senast visades eller uppdaterades.

**OBS!** Vid temperaturlarm växlar den visade temperaturen i användargränssnittet omedelbart till den faktiska lufttemperaturen inne i skakapparatens kammare för att skydda proverna på bästa möjliga sätt. Detta kan leda till ett hopp i den visade kammartemperaturen i diagramfunktionen.



Bild 73: Hastighets- och temperaturdiagram

Du kan spara data i hastighets- eller temperaturdiagram på ett USB-minne anslutet till skakapparatens USB-port genom att trycka på **Exportera diagramdata**.

### Redigera diagram

1. Tryck på **Redigera** för att redigera diagrammet.



Bild 74: Redigera diagram: Välj visning av temperatur eller hastighet

2. Tryck på **Temp** eller **Hastighet** för att välja vilka data som ska visas.

- 3. Tryck på nedpil för att visa alternativ för Datumintervall.
- 4. Tryck för att välja 1 dag, 7 dagar, eller Anpassa för valfritt datumintervall.
- 5. Dra värdena för år, månad och dag uppåt eller nedåt med fingret för att ställa in startdatum och tryck på **Ställ in startdatum**.
- Dra värdena för år, månad och dag uppåt eller nedåt med fingret för att ställa in slutdatum och tryck på Ställ in slutdatum.



Bild 75: Välja datumintervall för diagramvisning

- 7. Tryck på nedpil för att visa alternativ för Tidsintervall.
- 8. Tryck för att välja Hela dygnet, eller tryck på Anpassa för att välja valfri tidsperiod:
  - a. Dra värdena för timmar och minuter uppåt eller nedåt med fingret för att ställa in daglig starttid och tryck på **Ställ in starttid**.
  - b. Dra värdena för timmar och minuter uppåt eller nedåt med fingret för att ställa in daglig sluttid och tryck på **Ställ in sluttid**.

**OBS** Genom att välja start- och sluttid visas diagramdata för exakt denna tidsperiod. Om du har valt alternativet Hela dygnet börjar diagramdata registreras kl 00.00 och slutar kl 23.59 varje dag.



Bild 76: Välja tidsintervall för diagramdata

9. Tryck på **Spara** för att spara ändringar eller inställningar.

# 3.3. Tillbehör



Fysisk och biologisk skada orsakad av trasiga kärl. Felaktigt monterade tillbehör kan leda till krossat glas och utspillda prover.

SE UPP

Se till att tillbehören är korrekt monterade genom att använda rätt verktyg och skruvar.

Se till att tillbehören passar på plattformen.

Använd alltid kärl med tillbehör av passande storlek.



SE UPP

Skärskador från vassa kanter.

Var försiktig när du hanterar med plattformar och andra tillbehör.



SE UPP

Fingrar kan klämmas av plattformar i rörelse. Placera aldrig fingrarna på eller under en plattform medan den rör sig.

**OBS** Det är ditt ansvar att se till att tillbehören är korrekt monterade.

De monteringsprocedurer som beskrivs gäller för alla modeller av skakapparater som listas i denna bruksanvisning, tills annat anges.

Använd alltid lämpliga verktyg, i första hand verktygen som medföljer tillbehöret. Om du förlorar några av dessa delar kan du beställa nya reservdelssatser enligt lista ("Tillbehör" på sidan 17). Använd inte andra verktyg än de som föreskrivs, eller andra skruvar än de som medföljer.

# 3. 3. 1. Montera plattform

**SE UPP** En fullständig lista över kompatibla plattformar för varje skakapparat finns i avsnitt "1. 2. 1. Plattformar" på sidan 18. Temperaturstyrda skakapparater kan användas endast med den förinstallerade plattformen.

SE UPP Vassa kanter kan ge skärskador. När en plattform ska tas bort, ta tag i dess undersida. Lyft inte upp en plattform genom att hålla i kolvhållare monterade på den. Var försiktig när du hanterar med plattformar och andra tillbehör.

**SE UPP** Risk för klämskador från en rörlig 11 x 14-plattform med Solaris 4000 I/ 4000 R-modellerna. Använd inte 11 x 14-plattformen med Solaris 4000 I/4000 R.

SE UPP Använd endast de skruvar och verktyg som medföljer plattformen. Att använda andra skruvar eller fel verktyg kommer att leda till felaktig montering och kan därför medföra skador på skakapparat och tillbehör.

Montera alltid plattformar med **alla** skruvar.

Använd alltid rätt verktyg:

- 3/16"-verktyget (GT530066) för alla plattformar för Solaris 2000 och Solaris 2000 I/R och 4000 I/R
- 7/32"-verktyget (GT530080) för alla Solaris 4000-plattformars

## Universalplattformar

Solaris 2000 och 4000 levereras med en universalplattform, skruvar och verktyg. Vid behov kan ytterligare plattformar anskaffas. En fullständig lista över kompatibla plattformar för varje skakapparat finns i avsnitt "1. 2. 1. Plattformar" på sidan 18. 2000 I-, 2000 R-, 4000 I- och 4000 R-modellerna levereras med förinstallerad plattform. Varken plattform eller tillhörande skruvar medföljer som separata delar.



**Bild 78:** Universalplattform monterad på en Solaris 2000 (vänster) och en Solaris 4000 (höger) skakapparat

1. Placera försiktigt plattformen horisontellt över skakapparaten och dess infästningspunkter.

Plattformar för Solaris 2000 har 3 monteringshål. Plattformar för Solaris 4000 har 4 monteringshål.

Plattformar för Solaris 2000 I/R och 4000 I/R har tre monteringshål.

2. Dra åt skruvarna försiktigt för att fästa plattformen på skakapparaten. Sluta dra åt när verktyget börjar böjas. Tillbehör

### Dual Stack-plattformar

OBS Dual Stack-plattformar är inte kompatibla med Solaris 2000 I/R eller 4000 I/R.



Bild 79: Montera Dual Stack-plattform

Fäst den övre och den undre plattformen genom att montera dem med en pelare i varje hörn. Montera pelarna med rätt skruvar från den övre och den nedre plattformen.

Dra åt skruvarna försiktigt för att sätta ihop pelarna och plattformarna. Sluta dra åt när verktyget börjar böjas.

Den undre plattformen är den som har monteringshål för att ansluta den till skakapparaten. Se till att ha den undre plattformen nederst när du monterar.

#### Tillbehör

**OBS** Innan du placerar kärl på den monterade plattformen: Gör en slutlig kontroll för att vara säker på att plattformen inte vinglar.



Bild 80: Dual Stack-plattform på en Solaris 4000 skakapparat

1. Placera försiktigt den monterade Dual Stack-plattformen horisontellt över skakapparaten och dess monteringsplatta.

Plattformar för Solaris 2000 har 3 monteringshål, plattformar för Solaris 4000 har 4 monteringshål.

 Dra åt skruvarna försiktigt för att fästa plattformen på skakapparaten. Sluta dra åt när verktyget börjar böjas.

## 3. 3. 2. Montera kolvhållare och kärl

SE UPP Trasiga eller läckande kärl kan medföra biologisk skada. Felaktigt monterade tillbehör kan leda till att prover spills ut. Se till att tillbehören är korrekt monterade genom att använda rätt verktyg och skruvar. En PH2-skruvmejsel med skaftlängd 150 mm/6 in. eller längre krävs. (Artikelnummer 75004131). Se till att tillbehören passar på plattformen. Använd alltid kärl med tillbehör av passande storlek. Kärlen måste vara intakta och monterade på rätt sätt.

SE UPP Vassa kanter kan ge skärskador. Var försiktig när du hanterar med plattformar och andra tillbehör.

SE UPP Hållaren måste fästas med medföljande skruvar för att säker fixering ska garanteras. Montera alla medföljande skruvar i hållarens samtliga monteringshål.

**OBS** Om skruvarna är för långa riskerar plattformen att bli instabil, och dess funktion kan äventyras. Använd endast medföljande skruvar för fixeringen.

## <u>Kolvhållare</u>

Varje kolvhållare består av en hållare, en eller två fjädrar beroende på kolvhållarens/ kolvens storlek och skruvar för montering på plattformen. Använd endast de skruvar som medföljer kolvhållaren.



1) Fjäderrör 2) Fjäder 3) Skruvar 4) Ben 5) Fjäder

Bild 81: Detaljer, kolvhållare med 2 fjädrar

- 1. Om det behövs fäster du fjädern på benen som visas i illustrationerna.
- 2. Fjäderrören av gummi placeras mellan benen som visas på bilden. Vissa hållare använder två fjädrar. Den andra fjädern monteras runt basen på den hopsatta hållaren efter att den monterats på plattformen.
- 3. Fäst den hopsatta hållaren på plattformen med de medföljande skruvarna.

# <u>Kärl</u>

- Placera försiktigt det önskade kärlet i hållaren genom att först dra isär klämfjädern tillräckligt långt för att göra det möjligt att ställa in kärlet i hållaren. Skjut försiktigt in kärlet i rätt läge och fäst det i den bredare nederdelen av hållaren. Fjädern kommer att hålla kärlets hals säkert på plats.
- 2. Kontrollera att alla kärl sitter fast ordentligt innan du startar enheten.

När så är möjligt ska kärlet förseglas med en plugg för att förhindra att innehållet kastas ut vid blandningsprocessen.

## 3. 3. 3. Montera kvadratiska hållare

SE UPP Trasiga eller läckande kärl kan medföra biologisk skada. Felaktigt monterade tillbehör kan leda till att prover spills ut. Se till att tillbehören är korrekt monterade genom att använda rätt verktyg och skruvar. En PH2-skruvmejsel med skaftlängd 150 mm/6 in. eller längre krävs. (Artikelnummer 75004131). Se till att tillbehören passar på plattformen. Använd alltid kärl med tillbehör av passande storlek. Kärlen måste vara intakta och monterade på rätt sätt.

SE UPP Vassa kanter kan ge skärskador. Var försiktig när du hanterar med plattformar och andra tillbehör.



### Information om kvadratiska hållare

Varje kvadratisk hållare består av en bottenplatta med monteringshål och fyra fjäderbelastade ben. Dessutom medföljer skruvar för montering av hållaren på plattformen. Använd endast de skruvar som medföljer kolvhållaren.



1 Monteringshål 2 Fjäder 3 Ben 4 Skruvar

Bild 82: Information om kvadratiska hållare

- 1. Vid behov, haka fast alla fjädrar 2 på benen 3 (Bild 82).
- 2. Placera den hopsatta hållaren på plattformen och rikta in monteringshålen ①.
- 3. Fäst den hopsatta hållaren på plattformen med de medföljande skruvarna ④.
- 4. Sänk ned det kvadratiska kärlet ovanifrån.
- 5. Kontrollera att kärlet vilar på bottenplattan och är säkert förankrat i hållaren.

# 3. 3. 4. Montera provrörsställ

SE UPP Vassa kanter kan ge skärskador. Var försiktig när du hanterar med plattformar och andra tillbehör.

Varje provrörsställ består av två monteringskonsoler (metalltråd eller plåt) med en spärrknopp och två tumskruvar per provrörsställ för montering på plattformen. Varje provrörsställssats levereras monterad med ett provrörsställ och skumplastinlägg. Använd alltid de skruvar som medföljer stället.



Bild 83: Komplett provrörsställ med monteringskonsoler av metalltråd



- Provrörsställ
- ② Insatser av skumplast
- ③ Låsknopp
- ④ Monteringskonsol

Bild 84: Komplett provrörsställ med monteringskonsoler av plåt

- 1. Montera monteringskonsolerna på plattformen med tumskruvarna. Dra åt fingerskruvarna så långt det går med handkraft.
- 2. Sätt i provrörsstället i monteringskonsolerna.
- Ställ in önskad vinkel genom att använda låsknopparna. Dra åt låsknopparna så långt det går med handkraft.

# 3. 3. 5. Montera provrörsställhållare med justerbar vinkel

SE UPP Vassa kanter kan ge skärskador. Var försiktig när du hanterar med plattformar och andra tillbehör.

Provrörsställhållare och provrörsställ finns i olika storlekar (kapitel "1. 2. 5. Provrörsställhållare med justerbar vinkel" på sidan 28). Alla provrörsställhållare är justerbara i sju lägen från 0° vinkel i båda riktningarna, och kan spärras i 15°, 30° och 45°.

Använd alltid de skruvar som medföljer provrörsställhållaren.



Bild 85: Provrörsställhållare med justerbar vinkel med monterat provrörsställ

- 1. Montera provrörsställhållarna ① på skakapparatens plattform med medföljande skruvar.
- Dra metallflikarna ③ i den svängbara hållarens båda ändar något utåt ⑥ och sätt i provrörsstället ②.
- 3. Säkra provrörsstället 2 genom att försiktigt släppa flikarna 3.
- 4. Dra låsstiftens spärrknoppar (5) på hållarens båda gavlar utåt. Stiften frigörs genom vridning av knoppen ett fjärdedels varv.
- 5. Vrid provrörsställhållarens ① svängbara del ⑥ till önskad vinkel; 15°, 30° eller 45°.
- 6. Montera tillbaka låsstiftet (5) i positionshålet (4) och spärra stiftet genom att vrida knoppen ett fjärdedels varv.
- Ta bort provrörsstället genom att dra metallflikarna ③ i den svängbara hållarens båda ändar utåt ⑥ och lyft ut provrörsstället ①.

# 3. 3. 6. Montera hållare för mikrotiterplatta/deepwellplatta

SE UPP Vassa kanter kan ge skärskador. Var försiktig när du hanterar med plattformar och andra tillbehör.



- 1) Deepwellplatta
- ② Skruv
- ③ Hållare för
- mikrotiterplatta

Bild 86: Komplett mikrotiterplatta/deepwellplatta

- 1. Placera mikrotiterplattans ram på plattformen.
- 2. Montera mikrotiterplattans ram på plattformen med hjälp av skruvarna som medföljer mikrotiterplattan/deepwellplattan.
- 3. Sätt in mikrotiterplattan eller deepwellplattan i mikrotiterplattans ram.
- 4. Kontrollera att mikrotiterplattan eller deepwellplattan är säkert förankrad genom att försiktigt lyfta den.

# 3. 3. 7. Montera justerbar kärlhållare

SE UPP Vassa kanter kan ge skärskador. Var försiktig när du hanterar med plattformar och andra tillbehör.





① Monteringshål ② Justerbara stag ③ Låsrattar ④ Sidovägg ⑤ Glidsäker matta
 ⑥ Skruvar

Bild 87: Komplett justerbar kärlhållare

Tillbehör





① Monteringshål ② Justerbara stag ③ Låsrattar ④ Sidovägg ⑤ Glidsäker matta
 ⑥ Skruvar

Bild 88: Komplett kärlhållare justerbar i två höjder

Montering:

- 1. Lyft upp den glidsäkra mattan för att exponera monteringshålen (1) i Bild 88).
- 2. Placera den justerbara kärlhållaren på plattformen och rikta in monteringshålen.
- 3. Montera den justerbara kärlhållaren på plattformen med skruvarna 
  © som medföljer hållaren.
- 4. Placera den glidsäkra mattan 5 på hållarens botten.
- 5. Placera kärlet/kärlen på den glidsäkra mattan ⑤.
- 6. Lossa låsrattarna 3.
- 7. Justera stagen 2 så att kärlen är stadigt fixerade.
- 8. Dra åt låsrattarna 3.
- 9. Kontrollera att kärlet är stadigt fixerat genom att försiktigt lyfta det.

### 100 / 140

# 3. 3. 8. Montera bägarställ

SE UPP Vassa kanter kan ge skärskador. Var försiktig när du hanterar med plattformar och andra tillbehör.



Bild 89: Komplett bägarställ

- 1. Placera bottenplattan 2 på plattformen.
- 2. Placera bägarstället ① på bottenplattan.
- 3. Rikta in monteringshålen mot plattformen.
- 4. Fäst bägarstället med skruvarna ③ som medföljer produkten.

# 3. 3. 9. Montera separertratthållare

SE UPP Vassa kanter kan ge skärskador. Var försiktig när du hanterar med plattformar och andra tillbehör.

**OBS** Använd rätt skruvar. Två olika typer av skruvar medföljer. De längre skruvarna med röd gänglåsning används för det vertikala stativet. De kortare skruvarna utan gänglåsning används för montering av separertratthållaren på plattformen.

Separertratthållaren kan monteras på skakapparatens plattform på två sätt: horisontellt eller vertikalt.



Bild 90: Monteringslägen för separertratthållare

Det rakt horisontella läget ③ i Bild 90 kan justeras i höjdled till ett lutande läge; se ④. Separertratthållaren kan dessutom kombineras med ett vertikalt stativ (tillbehör) som möjliggör ett vertikalt läge; se ① och ②. Hållaren kan dessutom monteras med trattmynningen nedåt ①, eller omvänt för upp och nedvänd placering; se ②.

## Förbereda vertikal montering

I Bild 91 visas den kompletta separertratthållaren med ® vertikalt ® stativ.



Bild 91: Separertratthållare (vänster) med vertikalt stativ (höger)

- 1. Passa in separertratthållaren mot det vertikala stativets långa ände (1) och 2) i Bild 90).
- 2. Rikta in hållaren mot de fyra monteringshålen (2 i Bild 91) på det vertikala stativet:
  - a. Använd hålen för normal vertikal montering (③ i Bild 91) för montering med trattmynningen vänd nedåt.
  - b. Använd hålen för upp och nedvänd vertikal montering (④ i Bild 91) för montering med trattmynningen vänd uppåt.

## Montera på plattform

- 1. Placera den kompletta separertratthållaren på plattformen.
- 2. Rikta in monteringshålen mot plattformen.
- Fäst separertratthållaren på plattformen med skruvarna som medföljer hållaren (
   í Bild 91 och Bild 92).

# Montera tratten



Bakre tratthållare
Främre tratthållare
Låsratt för höjdjustering
Låsratt för längdjustering
Skruvar utan gänglåsning

Separertratt

Bild 92: Komplett separertratthållare

- 1. Lossa låsratten för längdjustering (5) och skjut den främre tratthållaren (3) till läget längst fram.
- 2. Sätt i trattens bakre ände ① i den bakre tratthållaren ②.
- Justera den främre tratthållaren ③ till bekväm höjd och fixera genom att dra åt låsratten för höjdjustering ④. Genom att modifiera höjdjusteringen kan separertratten lutas i önskat läge.
- 5. Fixera den främre tratthållaren ③ genom att dra åt låsratten för längdjustering ⑤.

# 3. 3. 10. Montera universalbricka

SE UPP Vassa kanter kan ge skärskador. Var försiktig när du hanterar med plattformar och andra tillbehör.



Bild 93: Universalbricka

Montering:

- 1. Placera universalbrickan 2 på plattformen.
- 2. Rikta in monteringshålen mot plattformen.
- 3. Fäst universalbrickan 2 med medföljande skruvar 3.
- 4. Placera gummimattan ① på brickan ②.
- 5. Placera flaskor på brickan och fixera med de medföljande O-ringarna ④ (Bild 94).

OBS O-ringar bör användas vid hastigheter över 100 v/min.



Bild 94: Universalbricka med ilagda behållare och O-ringar

**OBS** Var försiktig så att inte O-ringarna spänns ut för hårt när behållarna fixeras. När en O-ring spänns ut når den förr eller senare en punkt där den förlorar sin elasticitet och inte kan spännas ut mer. När denna punkt har uppnåtts, försök inte spänna ut O-ringen längre eftersom den då kommer att brista.

## 3. 3. 11. Montera och använda klistermatta

Klistermattor är lämpliga för behållare med bred och plan undersida, såsom Erlenmeyer-kolvar, Fernbach-kolvar, mediaflaskor, mikrotiterplattor, petriskålar, cellkulturkolvar, volymetriska kolvar och bägare. I allmänhet kan alla kärl som kan stå upprätt användas. Ju mindre kärlets basyta är i förhållande till dess höjd, desto lägre är emellertid maximal skakningshastighet. Höga behållare med liten bas är mindre lämpliga (Bild 95). Stora Erlenmeyer-kolvar (till exempel 3 000 ml) har starkare vidhäftning än små sådana (till exempel 100 ml).



Bild 95: Klistermatta: Lämpliga kolvformer

Två allmänna typer av klistermattor finns tillgängliga:

- Klistermatta med låg vidhäftning (transparent) 75004111 och 75004117
  - » enkel hantering tack vare lägre klibbighet
  - » särskilt lämpliga för bägare
  - » lämplig för alla kärl upp till 200 v/min (odling av vävnadskulturer)
- Klistermatta med hög vidhäftning (svart, transparent) 75004126 och 75004127
  - » tillförlitlig fixering av behållare tack vare högre klibbighet
  - » lämplig för alla kärl
  - » maximal hastighet 400 v/min (upp till 500 ml Erlenmeyer-kolvar)

Tillbehör

# <u>Verktyg</u>

Rollern som visas i Bild 96 är ett tillbehör för klistermattor. Den underlättar montering av klistermattan på plattformen.



Bild 96: Roller för klistermatta

### Montera torr klistermatta

- 1. Se till att skakapparatens plattform är ren och torr.
- 2. Dra bort skyddsfilmen från klistermattans ena sida.
- 3. Placera mattan med den vidhäftande sidan nedåt på plattformen.
- 4. Dra bort skyddsfilmen.

### Montera våt klistermatta

- 1. Dra bort skyddsfilmen från klistermattans båda sidor.
- 2. Blötlägg mattan i vatten eller skölj den noggrant.
- 3. Låt mattan rinna av.
- 4. Placera den fuktiga mattan på den rena plattformen.
- 5. Låt mattan torka på plattformen i 24 timmar före användning.

### Kärlegenskaper att beakta

Maximal hastighet är beroende av kärlets typ, dvs. storlek, form, basens form, material, ytjämnhet och renhet. Dessa egenskaper måste bedömas för att du ska kunna välja rätt klistermatta för dina behov.

I Tabell 53 framgår vilka material som är lämpliga för användning med klistermattor.

Plast/glas	Klibbighetsgrad	Metall	Klibbighetsgrad				
Glas (borsilikatglas)	5	Stål, rostfritt	5				
Polypropylen (PP)	3	Aluminium	5				
Sampolymerisad	3	Eloxerad	4				
polypropylen (PPCO)		aluminium					
Polyetylen (HPDE) (LPDE)	3						
Polykarbonat (PC)	4						
Polyetylentereftalat (PET)	4						
Glykolmodifierad	4						
polyetylentereftalat (PETG)							
Polymetylpenten (PMP)	3						
Polytetrafluoretylen (PTFE)	0*						
Sampolymerisad	0*						
tetrafluoretylen-							
hexafluorpropylen (FEP)							
Perfluoroalkoxypolymerer	2						
(PFA)							
Silikon	1*						
Gradering: 0 = ingen klibbighet, 5 = högsta klibbighet *Använd inte kärl graderade 0 eller 1							

(PTFE, FEP och silikon).

Tabell 53: Kompatibilitet för kärlmaterial

**OBS** Maximal skakhastighet kan uppnås endast med material som har klibbighetsgrad 5 i Tabell 53.

**OBS** För andra material än glas måste du göra egna test för att fastställa vilka material som kan användas med avseende på last, hastighet, körtid och temperatur.
#### Hastigheter och körtider för glaskärl (fyllt till 30 %)

Hastighets- och körtidsbegränsningar vid rumstemperatur för glaskärl (fyllt till 30 %). Förhöjd temperatur ger lägre klibbighet. Om temperaturen förändras kan kärlet lossna från klistermattan. Använd acklimatiserade kärl när skakningsprocessen startas.

	Mikrotiterplatta / Deepwellplatta	Bägare			
	96 x 2 ml	25 ml	125 ml	250 ml	500 ml
125 v/min	ingen körtidsbegränsn	iing			
150 v/min	72 timmar	72 timmar	72 timmar	72 timmar	72 timmar
200 v/min	72 timmar	72 timmar	72 timmar	72 timmar	72 timmar
250 v/min	72 timmar				
300 v/min					
350 v/min					
400 v/min					

Tabell 54: Klistermatta med låg vidhäftning – Hastighets- och körtidsbegränsningar vid rumstemperatur

	Mikroti- terplatta / Deepwell- platta	Erlenmeyer							
	96 x 2 ml	25 ml	125 ml	250 ml	500 ml	1000 ml	2000 ml	3000 ml	5000 ml
125 v/min	ingen körtids	sbegräns	ning						
150 v/min	72 timmar	72	72	72	72	72	72	72	72
		timmar	timmar	timmar	timmar	timmar	timmar	timmar	timmar
200 v/min	72 timmar	72	72	72	72	72	72	72	72
		timmar	timmar	timmar	timmar	timmar	timmar	timmar	timmar
250 v/min	72 timmar	72	72	72	72	72	72	48	24
		timmar	timmar	timmar	timmar	timmar	timmar	timmar	timmar
300 v/min	72 timmar	72	72	72	72	72	48	24	
		timmar	timmar	timmar	timmar	timmar	timmar	timmar	
350 v/min	48 timmar	48	48	48	48	48	24		
		timmar	timmar	timmar	timmar	timmar	timmar		
400 v/min	24 timmar	24	24	24	24				
		timmar	timmar	timmar	timmar				

Tabell 55: Klistermatta med hög vidhäftning – Hastighets- och körtidsbegränsningar vid rumstemperatur

#### Förberedelse

SE UPP Använd alltid skyddshandskar och skyddsglasögon vid hantering av glasprodukter. Kontrollera följande innan kärl placeras på klistermattan:

- 1. Kontrollera att kärl är oskadade; använd aldrig ett skadat kärl.
- 2. Rengör kärlet med en ren och torr duk så att all fukt avlägsnas.
- 3. Avlägsna allt fett och annan smuts.
- 4. Kontrollera att kärlets undersida och klistermattan är torra och rena.

SE UPP Även knappt synlig film eller en enda droppe vatten på kärlets undersida kan medföra svag vidhäftning och att kärlet lossnar från klistermattan efter en viss körtid.



Bild 97: Klistermatta: Torka kärlet torrt

5. Minska fyllvolymen eller försegla kärlet för att förhindra vätskespill.

SE UPP Även om ett kärl har perfekt vidhäftning kan en enda droppe vatten som rinner längs kärlets sida medföra svag vidhäftning och att kärlet lossnar från klistermattan efter en viss körtid.

#### Placera kärl

- 1. Placera kärlen på klistermattan och tryck fast dem.
- 2. Tryck eller dra kärlen försiktigt för att säkerställa tillförlitlig vidhäftning.

**OBS** Tänk på att kondensvatten kan bildas vid låga temperaturer eller långa körtider och medföra att kärlet lossnar från mattan. Prov- och kärltemperaturen får inte avvika från skakapparatens specificerade temperaturområde. Information om skakapparatens temperaturområde finns i "1. 1. Tekniska data" på sidan 11.

#### <u>Ta bort kärl</u>

- 1. Tryck på Stopp på skärmen eller vänta tills körningen är avslutad.
- 2. Vänta tills plattformen är helt stillastående.
- 3. Ta tag i kärlhalsen och dra kärlet åt sidan och vänta några sekunder. För kärl med stor bottenyta kan det ta några sekunder innan kärlet lossnar från klistermattan.



Bild 98: Klistermatta: Tryck eller dra kärlet försiktigt

SE UPP Tryck eller dra försiktigt; använd inte överdriven kraft! Kärlet kan lossna plötsligt så att innehållet spiller över!

4. Om ett kärl vägrar lossna kan du hälla lite vatten runt dess bas.

Låt klistermattan lufttorka innan den används på nytt.

#### Ta bort klistermattan

1. Ta bort klistermattan genom att lyfta upp ena sidan och långsamt dra loss mattan från plattformen (Bild 99).



Bild 99: Klistermatta: Lyfta och dra bort klistermattan

#### Rengöring och desinficering

Om klistermattans klibbighet minskar behöver mattan rengöras.

SE UPP Använd inte lösningsmedel eller rengöringsmedel annat än milt diskmedel utan hudvårdstillsatser eller liknande. Rekommenderad koncentration är högst två ml rengöringsmedel per liter vatten.

#### Tillbehör

**OBS** Om klistermattan kommer i direkt kontakt med koncentrerat rengöringsmedel minskar mattans klibbighet. I detta fall är det inte tillräckligt att endast skölja mattan. Skölj mattan, blötlägg den i vatten i 24 timmar och låt den sedan torka i 24 timmar för att den ska återfå full klibbighet.

OBS Desinficera klistermattan endast med 70 % isopropanol eller 70 % etanol.

#### Lätt nedsmutsad eller dammig matta:

- 1. Våttorka eller desinficera mattan på plattformen.
- 2. Fukta mattan.
- 3. Torka bort smutsresterna med en svamp.
- 4. Våttorka med en svamp genomdränkt med vatten.
- 5. Desinficera vid behov.
- 6. Låt torka i 24 timmar.

#### Kraftigt nedsmutsad matta:

- 1. Ta bort mattan från plattformen.
- 2. Blötlägg mattan i vatten med lite tillsatt diskmedel.
- 3. Gnugga mattan grundligt med en svamp för att avlägsna alla smutsrester.
- 4. Skölj mattan noggrant med vatten.
- 5. Desinficera vid behov.
- 6. Låt torka i 24 timmar.
- Montera mattan enligt beskrivningen i avsnitt "Montera torr klistermatta" på sidan 107 eller "Montera våt klistermatta" på sidan 107.

#### 3. 3. 12. Montera gasfördelare

Med gasfördelaren är det möjligt att spruta in inert gas såsom kväve eller koldioxid i flera kärl under skakningsprocessen (se kapitel "1. 2. 12. Allmänna tillbehör" på sidan 42 för beställningsinformation).

Gasfördelaren är ett förinstallerat tillval på Solaris 2000 I/R och 4000 I/R. Den är även tillgänglig för installation i efterhand av kund. Solaris 2000 I/R och 4000 I/R har en tillbehörsport på enhetens sida (kapitel "1. 4. Produktöversikt" på sidan 44) för matning av gas via labbets gasledning till fördelaren. Gasfördelaren har ett inlopp för gastillförsel och åtta utlopp till kärl.



Bild 100: Gasfördelare monterad i skakapparaten Solaris 4000 R

#### Montering:

- 1. Fäst gasfördelaren <sup>(2)</sup> på den bakre väggen på skakapparatens kammare med medföljande skruvar <sup>(5)</sup>.
- 2. Ta bort plastpluggen från tillbehörsportens genomföring (se kapitel "1. 4. Produktöversikt" på sidan 44).
- Gör ett korsformat snitt med en vass kniv i gummigenomföringens membran på kammarens insida.
- 4. Klipp till en bit flexibel slang i lämplig längd för användning som gasledning ①.
- 5. Anslut gasledningen ① till det räfflade inloppet ④ på gasfördelaren.
- 6. Kontrollera att gasledningen ① inte blockerar lockspärrarna, och led den genom tillbehörsportens genomföring till utsidan.

#### Tillbehör

- 7. Anslut gasledningen ① till labbets gaskälla.
- 8. Klipp till åtta slangbitar och fäst dem på utloppen ③ på gasfördelaren. Om fler än åtta kärl används kan "Y"-förgrenare användas; för färre än åtta kärl, blockera oanvända slangar så att gas inte släpps ut i onödan.
- 9. Använd lämpliga pluggar och tätningar för att koppla gasslangarna till kärlen.

## 3.4. Lastning och normal användning



Risk för brand på grund av utlösta kemiska reaktioner. Använd inte skakapparaten vid hastigheter som gör att innehållet i kärlen kastas ut

VARNING

Öka hastigheten långsamt. Prova med vatten innan du använder kemikalier.



## Säkerheten kan försämras genom felaktig lastning och skadade tillbehör.

SE UPP

- Se till att lasten (tillbehör och prover) är symmetriskt ordnad till mitten av plattformen. När du använder en dual stack-plattform, se till att den nedre plattformen bär mer belastning än den övre plattformen.
- Överbelasta aldrig skakapparaten. För maximal belastning se "Tekniska data" på sidan 11. Lasten består av vikten av plattformen, tillbehör och prover som monterats på skakapparaten.
- Kontrollera att tillbehören är korrekt monterade innan du använder skakapparaten. Följ instruktionerna i avsnitt "Tillbehör" på sidan 88.



SE UPP

#### Risk för klämskador orsakade av igenfallande lock.

 Skakapparatens lock kan oavsiktligt falla igen och klämma fingrar och händer. Detta riskerar att hända om lockfjädern är defekt och inte kan hålla locket i helt öppet läge. Använd inte skakapparaten med defekt lockfjäder. Kontakta Thermo Fisher Scientifics kundtjänst och beställ en ny lockfjäder och byt den omgående.

#### Lastning

Se till att belastningen, inklusive tillbehör och prover, arrangeras symmetriskt till mitten av plattformen. När du använder en dual stack-plattform och inte utnyttjar hela kapaciteten: för att uppnå bästa prestanda ska lasten främst placeras på den undre plattan och ordnas symmetriskt till mitten av plattformen.

Vid användning av temperaturstyrd skakapparat, se till att lasten placeras så att det finns tillräckligt med fritt utrymme när locket stängs, och att behållarna kan röra sig fritt med locket stängt så att kollision och efterföljande skada på behållarna undviks.



Bild 101: Exempel på en rätt laddad plattform

_												ē.		
•	٠	٠			*	٠	٠	*	٠	٠	- f		$\sim$	þ
•	٠	٠	٠	٠	٠	٠	*	*	٠	٠	. "	1940 1940	(•))¥	
•	٠	٠	٠	*	*	٠	*	*	*	*	*			
•		٠	*	٠	*	*	*	•	•			×		þ
•	٠	*	٠	*	*	*	*	۰	٠	*	٠	٠	·(·)·~	
•			*	٠	*	•		٠	•	*	•	٠		
•	٠	•	٠	,	•	*	٠	*	۰	*	٠	٠		þ
•		*				*	*	٠		*	*	e		
•	٠	•	٠	•	•	•	٠	*	٠	٠	٠	٠		
•				*		*		*	*	*	*	e		þ
Ŀ	•	•		•	*	٠	*	*	٠	•	٠	•		

Bild 102: Exempel på en dåligt laddad plattform

OBS Den maximala belastningen består av vikten av plattformen, tillbehör och prover som monterats på skakapparaten. Se "1. 1. Tekniska data" på sidan 11 för information om maximal last.

#### Normal användning

Information om begränsning av hastighet och belastning är endast en vägledning för produktanvändning. Smidig drift med låg eller ingen vibration är avhängig flera faktorer såsom fyllnadsgrad, typ och skick för kolvhållare och klistermattor, typ av kärl, placering av kärlen på plattformen, ytan som skakapparaten är placerad på och inställd hastighet.

När du startar en ny eller okänd kombination eller när någon av dessa parametrar ändras: Öka hastigheten över 175 rpm långsamt och kontrollera om det uppstår oönskade vibrationer i lasten eller om skakapparaten rör sig. Det är ditt ansvar att använda skakapparaten på ett säkert sätt.

Solaris-skakapparaterna är konstruerade för låg värmeavgivning och små vibrationer. Detta gör dem mycket lämpade för användning med många olika labbutrustningar såsom miljökammare, inkubatorer och labbkylar. På grund av att aktuella lokalförhållanden, inklusive hyllors och annan inrednings stabilitet, kan variera stort ansvarar användaren för säker drift vid användning med övrig labbutrustning. Öka hastigheten långsamt och var uppmärksam på oönskade instrumentrörelser.

#### Säker hastighet

Lastoberoende säker hastighet för Solaris-skakapparaterna framgår av Tabell 56. Lasten består av vikten av plattformen, tillbehör och prover som monterats på skakapparaten. Se "Tekniska data" på sidan 11 för information om maximal last.

Modell	Max. säker hastighet
Solaris 2000	175 v/min
Solaris 4000	175 v/min
Solaris 2000 I	250 v/min
Solaris 2000 R	250 v/min
Solaris 4000 I	250 v/min
Solaris 4000 R	250 v/min

Tabell 56: Maximal säker hastighet

#### Högsta hastighet

Med Solaris-skakapparaternas maximala hastighet 525 v/min är lastförmågan begränsad (Tabell 57). Lasten består av vikten av plattformen, tillbehör och prover som monterats på skakapparaten. Ta hänsyn till att hastighetsförmågan hos de kolvhållare och kärl som används kanske inte lämpar sig för den hastighet du vill använda.

Modell	Maximal säker last
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

Tabell 57: Maximal last med högsta hastighet

**OBS** Tänk alltid på att stigande hastighet gör att kolvhållarna börjar öppnas. Det medför tillfälliga vibrationer till systemet på grund av större rörelser i kärlen.

#### Vikt/hastighetskurvor

Nedanstående vikt/hastighetskurvor gör det lättare att ställa in en optimal kombination av hastighet och last vid normal användning. Det gröna området visar kombinationer av hastighet och belastning som inte orsakar några eller endast mycket små vibrationer vid skakning. Det röda området visar kombinationer som kan orsaka kraftiga vibrationer under skakningen och kan resultera i oönskad förflyttning av skakapparaten. Eftersom dessa diagram endast är till för vägledning, måste du vara uppmärksam när din kombination av hastighet och belastning kommer närmare det röda området.

Tänk på att dina specifika användningsförhållanden kan orsaka oönskat beteende hos last eller instrument innan du når den angivna gränsen. Öka hastigheten gradvis för att utforska beteendet hos din specifika belastning. Lasten består av vikten av plattformen, tillbehör och prover som monterats på skakapparaten.

OBS De blå linjerna i diagrammen avser tomvikt för en specifik plattform.

#### Solaris 2000



Bild 103: Solaris 2000 - normal användning



#### Solaris 4000



Lastning och normal användning



Bild 105: Solaris 2000 I / 2000 R - normal användning



#### Solaris 4000 I / 4000 R

Bild 106: Solaris 4000 I / 4000 R - normal användning

## 4. Underhåll och skötsel



VARNING

#### Risker vid hantering av farliga ämnen

- Vid skakning av farliga ämnen: Beakta
   Världshälsoorganisationen WHO:s handbok "Laboratory
   Biosafety Manual" och eventuella lokala bestämmelser.
   Vid skakning av mikrobiologiska prover från riskgrupp
   II (enligt Världshälsoorganisationen WHO:s handbok
   "Laboratory Biosafety Manual") måste aerosoltäta
   biologiska tätningar användas. Sök på internetsidan från
   Världshälsoorganisationen WHO (www.who.int) efter
   "Laboratory Biosafety Manual". För material i en högre
   riskgrupp måste extra säkerhetsåtgärder vidtas.
- Om toxiner eller patologiska substanser har kontaminerat skakapparaten eller dess delar måste lämpliga desinfektionsåtgärder vidtas ("Dekontaminering" på sidan 124, "Desinficering" på sidan 124).
- Om en farlig situation inträffar: Stäng av strömförsörjningen till skakapparaten och lämna området omedelbart.



VARNING

#### Hälsoskador från smittsamma ämnen

Om vätskor eller andra material spills ut och hamnar under plattformen: Stäng omedelbart av skakapparaten, dra ur kontakten och ta bort plattformen ("Plattformar" på sidan 18). Städa upp spillet enligt normala laboratorierutiner. Använd lämplig personlig skyddsutrustning.

Eventuella inre justeringar eller reparationer måste utföras av en auktoriserad servicetekniker. Skakapparatens kåpa får inte öppnas av användaren.

Följ instruktionerna i den produktinformation som medföljer tillbehöret. Där anges specifika detaljer om hur du underhåller och rengör det på rätt sätt. Använd följande information i det här kapitlet endast som riktlinje.

### 4.1. Allmän information

För personligt, miljömässigt och materiellt skydd, måste du rengöra och vid behov desinficera skakapparaten och dess tillbehör regelbundet.

Thermo Fisher Scientific rekommenderar att du rengör och desinficerar skakapparaten minst en gång i månaden. Normal inomhusluft innehåller tusentals cirkulerande mikroorganismer som kan bosätta sig i din skakapparat, och utsätta sätta dina odlingar för risker.

Thermo Fisher Scientific rekommenderar användning av 70% etanol, eller 70% isopropanol eller 10% eller mindre kvartärt ammoniumbaserat desinfektionsmedel.

**OBS** Mekanismen kan skadas om vätskor tränger in. Låt inte vätskor, särskilt inte några organiska lösningsmedel, komma i kontakt med mekanismen eller mekanismens lager. Organiska lösningsmedel bryter ner fettet i mekanismens lager. Om vätskor läcker in i delar av skakapparaten som inte är åtkomliga för rengöring, sluta använda skakapparaten och kontakta serviceavdelningen.

**OBS** lcke godkända förfaranden eller ämnen kan försämra material i skakapparat och leda till fel. Använd inga andra rengörings- eller dekontamineringsmetoder än de som beskrivs här, om du inte är säker på att de är lämpliga för de aktuella materialen. Använd endast rengöringsmedel som inte skadar utrustningen. Vid osäkerhet, kontakta tillverkaren av rengöringsmedlet.

OBS Med temperaturstyrda skakapparater ska locket lämnas öppet så att vätskor kan avdunsta.

- Dra ut nätsladden innan du rengör, desinficerar eller dekontaminerar.
- Ta bort monterade tillbehör och plattformar från skakapparat före rengöring, desinficering eller dekontaminering.
- Använd varmt vatten med ett milt rengöringsmedel med en mjuk trasa för att rengöra materialen. Om du är osäker: Kontakta Thermo Fisher Scientific. Skölj av med rent vatten och torka noggrant.
- Använd aldrig frätande rengöringsmedel såsom fosforsyra, blekmedel eller skurpulver.
- Använd endast desinfektionsmedel med ett pH-värde på 6-8.
- Rengör eventuellt spill omedelbart med en luddfri trasa, fuktad med ett icke frätande rengöringsmedel enligt tillverkarens anvisningar.
- Spill kan sippra in under plattformen. Om spill tränger in under plattformen ska denna tas bort och spillet på tråget under plattformen torkas upp. Desinficera eller dekontaminera vid behov.
- Kontrollera skakapparatens delar och ta bort utspillda tillväxtmedia eller skräp.

#### Inspektera tillbehör

OBS Använd inte någon skakapparat eller några tillbehör som visar tecken på skada. Vi rekommenderar att du regelbundet inspekterar tillbehören som en del av ditt rutinunderhåll för att garantera säkerheten.

Efter noggrann rengöring av tillbehören måste de inspekteras för skador, slitage och korrosion.

#### <u>Metalldelar</u>

I händelse av skada, såsom korrosion, slitage eller sprickor, får inte tillbehören användas mer.

#### <u>Plastdelar</u>

Kontrollera om det finns tecken på krackelering, blekning, blåmärken eller sprickbildning. Vid skada måste tillbehöret omedelbart tas ur drift.

#### **Klistermattor**

Kontrollera före varje användning att klistermattorna inte uppvisar tecken på slitage eller skador såsom krackelering, missfärgning, gulning, sprödhet, deformation, utnötning eller kemiska angrepp. Kassera omedelbart klistermattor som uppvisar tecken på förslitning eller skada.

#### 4.2. Rengöring

**OBS** Innan du använder en viss rengöringsmetod bör du kontrollera med tillverkaren av rengöringsmedlet att den föreslagna metoden inte kommer att skada utrustningen.

OBS Mekanismen kan skadas om vätskor tränger in. Låt inte vätskor, särskilt inte organiska lösningsmedel, komma på mekanismen eller mekanismens lager. Organiska lösningsmedel bryter ner fettet i mekanismens lager. Rengör enhetens utsida med en mjuk trasa indränkt med en mild tvål- och vattenlösning, skölj med rent vatten och torka torrt.

Se "Allmän information" på sidan 122 för rengöring av skakapparat och tillbehör.

OBS Med temperaturstyrda skakapparater ska locket lämnas öppet så att vätskor kan avdunsta.

#### Pekskärm

Rengöra pekskärmen:

- 1. Dra ur stickproppen.
- 2. Rengör pekskärmen med en torr mikrofiberduk.
- 3. Vid behov, fukta mikrofiberduken med vatten och torka av pekskärmen igen.
- 4. Låt ytorna lufttorka

## 4.3. Desinficering

VARNING Risker vid hantering av farliga ämnen. Vidrör inte infekterade delar. Farlig smitta är möjlig vid beröring av förorenade delar. Smittsamt material kan komma in i skakapparaten när ett kärl går sönder eller som ett resultat av spill. Se till att ingen människa utsätts för risker i händelse av kontamination. Desinficera omedelbart utsatta delar.

**OBS** Utrustningen kan skadas av olämpliga desinfektionsmetoder eller medel. Försäkra dig om att desinfektionsmedlet eller desinfektionsmetoden inte ger upphov till skador på utrustningen. Om du är osäker, kontakta tillverkaren av rengöringsmedlet. Följ säkerhetsoch användningsanvisningarna för det aktuella desinfektionsmedlet.

- 1. Torka alla delar och områden med 70% etanol enligt krav för den nivå av desinfektion du behöver. Fukta inte områden med exponerad elektronik.
- 2. Låt ytorna lufttorka.

OBS Med temperaturstyrda skakapparater ska locket lämnas öppet så att vätskor kan avdunsta.

Du är ansvarig för att desinficering utförs enligt de krav som gäller på din arbetsplats.

### 4.4. Dekontaminering

VARNING Risker vid hantering av farliga ämnen. Vidrör inte kontaminerade delar. Exponering för kontaminering är möjlig när man vidrör de förorenade delarna. Förorenat material kan komma in i skakapparaten när ett rör bryts eller som ett resultat av spill. Se till att ingen människa utsätts för risker i händelse av kontamination. Dekontaminera genast alla kontaminerade delar.

**OBS** Utrustningen kan skadas av olämpliga dekontamineringsmetoder eller medel. Försäkra dig om att dekontamineringsmedlet eller dekontamineringsmetoden inte ger upphov till skador på utrustningen. Kontakta tillverkaren av dekontamineringsmedlet om du är osäker. Följ säkerhets- och användningsanvisningarna för det aktuella dekontamineringsmedlet.

Följande metod föreslås av Thermo Fisher Scientific.

- 1. Torka alla delar och ytor med 70% etanol. Fukta inte områden med exponerad elektronik.
- 2. Låt ytorna lufttorka

OBS Med temperaturstyrda skakapparater ska locket lämnas öppet så att vätskor kan avdunsta.

Du är ansvarig för att dekontaminering utförs enligt de krav som gäller på din arbetsplats.

## 4.5. Autoklavering

**OBS** Överskrid aldrig de tillåtna värdena för autoklaveringstemperatur och -tid. Kemiska tillsatser i ångan är inte tillåtna.

Före autoklavering:

- Demontera alla tillbehör.
- Skölj noggrant bort alla spår av kemikalier eller rengöringsmedel. Avsluta med att skölja med destillerat vatten.

Följ denna tabell för att kontrollera om det går att autoklavera delarna:

Artikel	Autoklave- ringsbar	Specifikationer
Skakapparat	Nej	-
Plattform	Ja	121 °C, 20 min
Kolvhållare (utan fjäderrör)	Ja	121 °C, 20 min
Skruvmejsel	Nej	-
Hållare för mikrotiterplatta/deepwellplatta	Ja	121 °C, 20 min
Klistermatta	Nej	-
Bägarställ	Ja	121 °C, 20 min
Separertratthållare	Ja	121 °C, 20 min
Provrörsställ – Metalldelar	Ja	121 °C, 20 min
Provrörsställ – Plastdelar	Nej	-
Provrörsställ – Skumplastinlägg	Nej	-
Justerbar kärlhållare – Metalldelar	Ja	121 °C, 20 min
Justerbar kärlhållare – Glidsäker matta	Nej	-
Universalbricka – Metallbricka	Ja	121 °C, 20 min
Universalbricka – Gummimatta	Nej	-
Provrörsställhållare med justerbar vinkel – Metalldelar	Ja	121 °C, 20 min
Provrörsställhållare med justerbar vinkel – Plastdelar	Nej	-
Provrörsställhållare med justerbar vinkel – Skumplastinlägg	Nej	-

 Tabell 58: Möjlighet att autoklavera material

Säkerställ nödvändig sterilitet i enlighet med aktuella krav.

Utseende och färg kan ändras något efter autoklavering.

## 4.6. Temperaturkalibrering

Temperaturbörvärdet hos en temperaturstyrd skakapparat kan kalibreras med en mätutrustning som uppvisar hög precision. Därmed kan fullständig reproducerbarhet i den inkuberade (uppvärmda) eller kylda (kylda och uppvärmda) skakningsprocessen säkerställas.

#### 4. 6. 1. Etablerad praxis för temperaturkalibrering

Följ instruktionerna noggrant före och under temperaturkalibreringen. Vid störning eller avvikelse från rekommenderad praxis riskeras att provtemperaturen inte stabiliseras och kalibreringen misslyckas.

Etablerad praxis inkluderar:

- Använd en kolv med 250 ml volym.
- Fyll kolven med exakt 100 ml vatten.
- Stäng locket helt och håll det stängt under hela kalibreringsprocessen.
- Avvakta tillräckligt länge för att provtemperaturen ska hinna stabiliseras helt.
- Se till att omgivningsförhållanden såsom temperatur och luftfuktighet inte ändras under kalibreringsprocessen. Till och med små förändringar som att öppna en dörr eller ett fönster eller ändring av luftkonditioningssystemets inställningar kan äventyra kalibreringsresultatet.
- Vi rekommenderar att skakapparaten kalibreras årligen.

#### 4. 6. 2. Procedur för temperaturkalibrering

Utför kalibreringen så här:

- 1. Kalibreringen kan startas på tre olika sätt:
  - » Efter att du har slutfört den initiala startprocessen som beskrivs i "Första start" på sidan 54 ges möjlighet att utföra temperaturkalibrering istället för att avsluta startprocessen.
  - » Skakapparaten visar en påminnelse om årlig kalibrering.
  - » Du kan starta kalibreringen genom att trycka på Inställningar i navigeringsfältet. Tryck sedan på Kalibrering på skärmen Temp. kalibrering.
  - » Du kan vid behov trycka på Inställningar i navigeringsfältet och sedan på Kalibrering för att starta processen.

OBS Du kan inaktivera kalibreringspåminnelser genom att markera kryssrutan **Inaktivera** kalibreringspåminnelser på skärmen **Inställningar**.

2. Vänta tills skärmen Temp. kalibrering (Bild 107) visas.



Bild 107: Kalibrering: Startskärmen

3. Tryck på Starta kalibrering.



Bild 108: Kalibrering: Uppmaning att fylla på Erlenmeyer-kolv

- 4. Följ instruktionerna på skärmen: Fyll 250 ml Erlenmeyer-kolven med 100 ml vatten.
- 5. Placera kolven i en hållare och fixera den ungefär i mitten av skakapparatens plattform.
- 6. Sänk ned sonden för en mätutrustning som har hög precision i kolven. Sonden ska vara nedsänkt i vätskan utan att vidröra kolvens botten eller sidor.
- 7. Försegla kolven så att ingen vätska spills ut.
- 8. Stäng skakapparatens lock.

9. Tryck på Nästa för att fortsätta.



Bild 109: Kalibrering: Uppmaning att starta kalibreringen

OBS Skakningsprocessen startar omedelbart när du trycker på Start.

 Följ instruktionerna på skärmen: Tryck på Start för att använda den förinställda temperaturen, eller tryck på Temperaturbörvärde och ange önskad temperatur innan du trycker på Start.

På skärmen visas hur skakapparatens kammare värms upp (eller kyls ned) till den valda kalibreringstemperaturen.



Bild 110: Kalibrering: Kammaren stabiliseras till kalibreringstemperatur

11. Avvakta minst 150 minuter för att kammartemperaturen ska uppnå ett stabilt tillstånd.

Temperaturen är stabiliserad när meddelandet **Kammartemperaturen är** stabiliserad (Bild 111).



Bild 111: Kalibrering: Kammaren har stabiliserats till kalibreringstemperatur

- 12. Följ instruktionerna på skärmen:
  - » Så snart skärmen i Bild 111 visas, jämför upprepade gånger temperaturavläsningen för termoelementet. När du bedömer att temperaturen är stabil kan du antingen trycka på **Redigera** för att korrigera värdet, eller på **Avsluta** för att slutföra kalibreringsprocessen.



Bild 112: Kalibrering: Korrekt temperatur visad av skakapparaten

- 13. Om temperaturen behöver korrigeras kan du dra värdet som visas på skärmen uppåt eller nedåt med fingret tills det överensstämmer med den temperatur som har uppmätts med mätutrustningen.
- 14. Tryck på Spara för att spara ändringen.

15. Meddelandet Temperaturkalibrering klar bekräftar att processen är slutförd.

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

Bild 113: Kalibrering: Slutförd temperaturkalibrering

Vid behov kan du trycka på Kalibrera om för att göra om kalibreringen.

#### 4. 6. 3. Temperaturkalibreringsvärden

Du kan när som helst visa den aktuella kalibreringstemperaturen:

- 1. Tryck på **Inställningar** i navigeringsfältet. Tryck sedan på **Kalibrering** på skärmen **Temp. kalibrering**.
- Skärmen Temp. kalibrering i Bild 107 visas. På denna skärm visas aktuell Kalibreringstemperatur.

OBS Med temperaturstyrda skakapparater kan du visa kalibreringstemperaturen även på skärmen för temperaturinställning. Ett exempel visas i avsnittet "Ställa in temperatur" på sidan 64.

#### 4.7. Installera inbyggd programvara

När du trycker på Inbyggd programvara på skärmen Filer och information ("Filer och information" på sidan 77) måste du ange administratörslösenordet för att visa skärmen Inbyggd programvara (Bild 114). Härifrån kan du installera uppdaterad inbyggd programvara från ett USB-minne.

Beroende på ditt serviceavtal kan en auktoriserad servicetekniker tillhandahålla en länk till en ny version av den inbyggda programvaran. Installera inbyggd programvara endast från en auktoriserad källa. Programvara från oauktoriserade källor kan ogiltiggöra Thermo Fisher Scientifics garanti.

- 1. Hämta zip-filen från länken som har erhållits från serviceteknikern.
- 2. Extrahera zip-filen till ett tomt USB-minne.

OBS Kontrollera att USB-minnet har tillräckligt med utrymme för extrahering av zip-filen.

- 3. Om skakapparaten är igång, stoppa den genom att trycka på **Stopp** på skärmen.
- 4. Kvittera eventuella larm; se beskrivningen i avsnittet "Status" på sidan 66.

OBS Om du försöker uppdatera programvaran medan en skakningsprocess pågår visas meddelandet **Uppdatering av skakapparaten kan inte göras medan en skakningsprocess pågår.** Tryck på knappen **Stoppa skakapparaten** i meddelandet för att fortsätta uppdateringen.

- 5. Klicka på ikonen Filer och information på navigeringsfältet.
- 6. Rulla ned till slutet av skärmen **Filer och information** tills knappen **Inbyggd programvara** visas.
- 7. Tryck på **Inbyggd programvara** för att visa skärmen Inbyggd programvara med nuvarande programvaruversion.
- 8. Sätt i USB-minnet i en av skakapparatens USB-portar. USB-portarnas placering beskrivs i avsnittet "Produktöversikt" på sidan 44.
- 9. Ange administratörslösenordet.



Bild 114: Skärmen Inbyggd programvara

10. Tryck på Uppdatera inbyggd programvara.

**OBS** Om inget USB-minne hittas visas meddelandet **Sätt i ett USB-minne med den inbyggda programvara som du vill installera för att starta uppdateringen.** Sätt i USB-minnet, alternativt kontrollera om det är korrekt anslutet till USB-porten. 11. Om det finns någon ny version av den inbyggda programvaran visas denna på skärmen Uppdatera inbyggd programvara (Bild 115). Tryck på den inbyggda programvara som du vill uppdatera.

Firmware Upgrade	×	Firmware Upgrade	×
Select firmware to upgrade		Select firmware to up	ograde:
O H.M.I.		Он.м.	
Main Controller		Main Controller	
O Parameters		O Parameters	
Select		Select	

Bild 115: Välja inbyggd programvara att installera

OBS Inbyggd programvara som är aktuell är gråtonad och kan inte väljas. Om det inte finns någon inbyggd programvara alls att uppdatera visas meddelandet **Inga** programuppdateringar finns för tillfället.

- 12. Tryck på Välj.
- Om USB-minnet innehåller mer än en version av den inbyggda programvaran uppmanas du att välja en version. Tryck på önskad programversion och tryck sedan på Välj.



Bild 116: Välja version av inbyggd programvara för installation

**OBS** Om du är osäker på vilken version du ska installera, kontakta serviceteknikern som har tillhandahållit länken till programvaran.

14. De programversioner som kommer att uppdateras visas på skärmen. Tryck på **OK** för att starta uppdateringen.



Bild 117: Lista med programversioner som ska uppdateras

**OBS** Installationsprocessen är automatisk; när du väljer en inbyggd programvara för uppdatering så hämtar installationsrutinen automatiskt alla tillhörande programkomponenter som krävs för uppdateringen.

**OBS** Om du väljer alternativet HMI i listan (Bild 115) uppdateras endast skakapparatens grafiska användargränssnitt. Pekskärmen kommer då att vara inaktiverad i upp till en minut efter påbörjad uppdatering. Stäng <u>inte</u> av skakapparaten under denna tid; i annat fall slutförs inte uppdateringen korrekt. Efter en kort stund startar skakapparaten om och är därefter klar att användas.

**OBS** Om du väljer två eller fler komponenter i listan (Bild 115) för uppgradering av <u>flera</u> programkomponenter samtidigt får skakapparaten <u>inte</u> stängas av medan uppdateringen pågår.

15. Olika statusmeddelanden om programuppdateringen visas under installationen (Bild 118).



Bild 118: Statusmeddelanden under installation av uppdaterad inbyggd programvara

16. När du uppmanas att starta om skakapparaten, stäng av strömbrytaren och slå sedan på den igen.

## 4.8. Plattformsbyte

Den grundläggande proceduren för montering av plattform beskrivs i kapitel "3. 3. 1. Montera plattform" på sidan 89. En fullständig lista över extra plattformar för varje skakapparat finns i avsnitt "1. 2. 1. Plattformar" på sidan 18.

**OBS** Temperaturstyrda skakapparater kan användas endast med den förinstallerade plattformen.

#### 4.9. Service

Thermo Fisher Scientific rekommenderar att skakapparaten och tillbehören servas en gång per år av en auktoriserad servicetekniker. Kundtjänstteknikern kontrollerar följande:

- den elektriska utrustningen
- uppställningsplatsens lämplighet
- säkerhetssystem
- använda tillbehör
- fastsättning av hållare och plattformar och andra tillbehör på skakapparaten
   Före service ska skakapparat och tillbehör noggrant rengöras och dekontamineras för att säkerställa att fullständig och säker inspektion kan utföras.

För dessa funktioner erbjuder Thermo Fisher Scientific inspektions- och serviceavtal. Nödvändiga reparationer utförs kostnadsfritt inom ramen för garantivillkoren, och därutöver mot betalning. Detta gäller endast om skakapparaten har servats av en servicetekniker som auktoriserats av Thermo Fisher Scientific.

### 4.10. Leverans och avyttring

VARNING Hälsoskador från smittsamma ämnen. När skakapparaten och tillbehören tas ur drift måste du rengöra och vid behov desinficera eller dekontaminera dem. Vid tvivel, kontakta Thermo Fisher Scientific kundtjänst.

För avfallshantering av skakapparaten gäller reglerna i ditt land. Kontakta Thermo Fisher Scientific kundtjänst för avfallshantering av skakapparaten. Kontaktinformation finns på sista sidan i denna manual. Du kan också besöka <u>www.thermofisher.com</u>.

Följ instruktionerna om transport och frakt ("Transportera" på sidan 52, "Leverans" på sidan 59).

## 5. Felsökning

Nr.	Beskrivning	Lösningar	lkon
Fel			
11999	Ett internt fel har inträffat	Ett internt fel har inträffat. Starta om skakapparaten genom att stänga av strömbrytaren och sedan slå på den igen. Om felmeddelandet fortfarande visas, kontakta en servicetekniker.	
Meddeland	len		
3	Temperaturkalibrering behövs. Den senaste gjordes för ett år sedan.	Ett år har passerat sedan temperaturkalibrering senast gjordes. Utför temperaturkalibrering enligt avsnitt "4. 6. Temperaturkalibrering" på sidan 126.	CAL
4	Strömmen återupptas och automatisk omstart aktiverad.	Ett strömavbrott inträffade under den senaste körningen. När strömmen kom tillbaka återupptogs körningen automatiskt.	Ŧ
6	Givare för omgivningstemperatur ur funktion.	Temperaturen som är uppmätt av omgivningstemperaturgivaren är osannolik.	
10	Hastigheten för fläkt 1 överstiger gränsvärdet.	Hastigheten för fläkt 1 är orimligt hög.	•2
11	Hastigheten för fläkt 2 överstiger gränsvärdet.	Hastigheten för fläkt 2 är orimligt hög.	•2
12	Hastigheten för fläkt 3 överstiger gränsvärdet.	Hastigheten för fläkt 3 är orimligt hög.	•2
13	Hastigheten för fläkt 4 överstiger gränsvärdet.	Hastigheten för fläkt 4 är orimligt hög.	•2
20	Locket har varit öppet under lång tid.	Stäng skakapparatens lock så att kammartemperaturen bibehålls.	
Larm			
1	Larm om hög temperatur	Kammartemperaturen överskrider övre gränsvärde. Kontrollera prover och omgivningsförhållanden och/eller ändra aktuella inställningar.	
2	Larm om låg temperatur	Kammartemperaturen underskrider undre gränsvärde. Kontrollera prover och omgivningsförhållanden och/eller ändra aktuella inställningar.	H
3	Automatisk omstart efter strömavbrott misslyckades.	Ett strömavbrott inträffade under den senaste körningen. Det gick inte att utföra automatisk omstart.	Ç,
5	Accelerationen är för långsam. Önskad inställd hastighet kan inte nås.	Den önskade inställda hastigheten kunde inte nås inom tidsgränsen. Kontrollera dina inställningar och/eller belastning (tillbehör och prover) på plattformen.	$\odot$

#### Leverans och avyttring

Nr.	Beskrivning	Lösningar	Ikon
7	Hastighetsmätning för fläkt 1 rapporterar oväntat stillestånd.	Stäng av enhetens strömbrytare och slå sedan på den igen. Om felmeddelandet fortfarande visas, kontakta en servicetekniker.	•2
8	Hastighetsmätning för fläkt 2 rapporterar oväntat stillestånd.	Stäng av enhetens strömbrytare och slå sedan på den igen. Om felmeddelandet fortfarande visas, kontakta en servicetekniker.	•2
9	Hastighetsmätning för fläkt 3 rapporterar oväntat stillestånd.	Stäng av enhetens strömbrytare och slå sedan på den igen. Om felmeddelandet fortfarande visas, kontakta en servicetekniker.	•2
10	Hastighetsmätning för fläkt 4 rapporterar oväntat stillestånd.	Stäng av enhetens strömbrytare och slå sedan på den igen. Om felmeddelandet fortfarande visas, kontakta en servicetekniker.	•2
12	Startfel - ingen hastighetssignal tas emot.	Skakapparatens plattform är blockerad. Kontrollera om det finns tillräckligt med utrymme runt plattformen och/eller minska belastningen (tillbehör och prover) på plattformen. Tryck därefter på START igen på skakapparaten. Om felmeddelandet fortfarande visas, kontakta en servicetekniker.	
24	Hastighetsmätningsfel under körning.	Onormal hastighetsändring har upptäckts. Kontrollera belastningen (tillbehör och prover) och/eller kolvhållare på plattformen. Tryck därefter på START igen på skakapparaten. Om felmeddelandet fortfarande visas, kontakta en servicetekniker.	-
26	Hastighetsmätningen upptäckte oväntat stillestånd under kömingen.	Se till att det finns tillräckligt med utrymme runt plattformen och/eller minska belastningen (tillbehör och prover) på plattformen. Kontrollera om säkringen på baksidan av skakapparaten har löst ut ("Säkring" på sidan 47). Tryck därefter på START igen på skakapparaten. Om felmeddelandet fortfarande visas, kontakta en servicetekniker.	$\bigcirc$
82	Motorströmmätning detekterar överbelastning	Överström har upptäckts på motorn. Förbjudet att lasta eller tömma plattformen under körning. Sänk hastigheten eller justera belastningen (tillbehör och prover) på plattformen.	<b></b> 0
83	Motorströmmätning utanför gränserna	Överström har upptäckts på motorn. Förbjudet att lasta eller tömma plattformen under körning. Sänk hastigheten eller justera belastningen (tillbehör och prover) på plattformen.	0

Tabell 59: Lista över fel, felmeddelanden och larm

**OBS** Om ett felmeddelande visas som inte finns med i denna tabell, kontakta en servicetekniker.

## GPL (General Public License)

Delar av enhetens programvara använder öppen källkod publicerad under GPL, LGPL eller annan licens för öppen källkod. Dessa bibliotek beskrivs i tabellen nedan. De aktuella bibliotekens källkod (från tredje part) kan erhållas från Thermo Fisher Scientific i de fall den relevanta licensen medger detta. Licensvillkoren för den aktuella öppna källkoden ingår i paketet med den öppna källkoden.

#### Bibliotek från tredje part

Bibliotek	Version	Spridning via	Licens
Qt	5.8	BSP	LGPLv3
Log4Cplus	1.2.0	Programvara	ApachePublicLicensev2/förenkladBSD-typlicens
boost	1.72.0	Programvara	Boost-licens 1.0
json11	1.0.0	Programvara	MIT-licens

\_\_\_\_\_

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

## Index

#### <u>A</u>

Anpassningsmenyn 76 Anslutningar 46 Autoklavering 125 Automatisk omstart 73 Avyttring 134

#### В

Behörighetskontroll 71

#### D

Datum 74 Datumformat 75 Dekontaminering 124 Desinficering 124 Diagram 86 Direktiv 43 Display 73 Drift 60 Dual Stack-plattformar 91

#### Ε

Enhetsnamn 76 Ethernet 47

#### F

Felsökning 135 Filer och information 77

#### G

GPL (General Public License) 137 Grafiskt användargränssnitt 60

#### Η

Hastighetsinställningar 73 Händelselogg 84 Högsta hastighet 118

Inspektera tillbehör 123 Inställningar 68

#### Κ

Kalibrering 72 Kontroller 73 Korrekt användning 5

#### L

Larm 69 Lastning 115, 116 Leverans 59, 134 Leveransomfattning 49 Ljusstyrka 73

#### Μ

Meddelanden 69 Metalldelar 123 Montera bägarställ 101 Montera gasfördelare 113 Montera hållare för mikrotiterplatta/ deepwellplatta 98 Montera justerbar kärlhållare 99 Montera kolvhållare och kärl 92 Montera kvadratiska hållare 94 Montera och använda klistermatta 106 Montera plattform 89 Montera provrörsställ 96 Montera provrörsställhållare med iusterbar vinkel 97 Montera separertratthållare 102 Montera universalbricka 105

#### Ν

Normal användning 115, 117 Nätanslutning 47

#### Ρ

Plastdelar 123

Plattformar 18 Produktöversikt 44 Program 78

#### R

Region 76 Rengöring 123

#### S

Service 77, 134 Signalord och symboler 6 Skötsel 121 Snooze Timeout 70 Specifikationer 11 Språk 74 Standarder 43 Start 54 Ställa in temperatur 64 Symboler som används i bruksanvisningen 7 Säker hastighet 117 Säkring 47

#### Т

Tekniska data 11 Temperaturbörvärde 64, 73 Temperaturkalibrering 126 Tid 75 Tidsläge 73 Tillbehör 17, 88 Transportera 52

#### U

Underhåll 121 Universalplattformar 90 Uppställningsplats 50 USB 48

#### V

Viloläge 75 Volym 69

# thermo scientific





Tillverkad för

Thermo Elektron LED GmbH Zweigniederlassung Osterode Am Kalkberg, 37520 Osterode am Harz Germany Ursprungsland: 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 är den ursprungliga bruksanvisningen. Denna bruksanvisning är en översättning av den ursprungliga bruksanvisningen.

#### thermofisher.com/shaker

© 2019–2024 Thermo Fisher Scientific Inc. Eftertryck och all annan kopiering förbjuden.

Alla varumärken tillhör Thermo Fisher Scientific Inc. och dess dotterbolag om inte något annat anges. Inte alla produkter finns tillgängliga i alla länder. Du kan få närmare information på förfrågan från din lokala försäljningspartner.

Bilderna som finns publicerade i denna bruksanvisning är bara till för referens. Där visade inställningar och språk kan avvika.

Australien +61 39757 4300

Österrike +43 1 801 40 0

Belgien +32 53 73 42 41

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

Frankrike +33 2 2803 2180

**Tyskland nationellt, avgiftsfritt** 0800 1 536 376

**Tyskland international** +49 6184 90 6000 Indien +91 22 6716 2200

**Italien** +39 02 95059 552

**Japan** +81 3 5826 1616

Nederländerna +31 76 579 55 55

**Nya Zeeland** +64 9 980 6700

Nordeuropa/Baltikum/OSS +358 10 329 2200

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

Andra asiatiska stater +852 2885 4613

**Andra länder** +49 6184 90 6000

