

Thermo Scientific Solaris sorozatú rázógép

Használati utasítások

70900205-d • 2023-03



Tartalom

Előszó		5
A kézikön	yvről	. 5
Rendeltet	ésszerű használat	. 5
	Nyitott rázógép	5
	Inkubált rázógép (csak fűtött) és hűtött rázógép (hűtött és fűtött)	5
	Szükséges tudnivalók	5
Jelek és k	ifejezések	. 6
	A rázón és tartozékain elhelyezett jelek	6
	A kezelési útmutatóban használt jelek	7
Biztonság	i előírások	. 7
1. Műszaki	adatok	11
1. 1. Műs	zaki adatok	11
	Thermo Scientific Solaris 2000 rázó	. 11
	Thermo Scientific Solaris 4000 rázó	. 12
	Thermo Scientific Solaris 2000 I rázó	. 13
	Thermo Scientific Solaris 4000 I rázó	. 14
	Thermo Scientific Solaris 2000 R rázó	. 15
	Thermo Scientific Solaris 4000 R rázó	. 16
1. 2. Tart	ozékok	17
1. 2. 1.	Tálcák	18
1. 2. 2.	Lombiktartó	19
1. 2. 3.	Lombiktartó Mikrotiter/Deepwell tálcákhoz	23
1. 2. 4.	Kémcsőállványok	24
1. 2. 5.	Állítható szögű kémcsőállványok	28
1. 2. 6.	Állítható edénytartó	30
1. 2. 7.	Szögletes edénytartó	32
1. 2. 8.	Nalgene főzőpohár állványok	34
1. 2. 9.	Szétválasztó tölcsértartó	38
1. 2. 10		39
1. 2. 11	Untapados szonyegek	40
1. 2. 12		42
1.3. Sza	bvanyok es irányelvek	43

1. 4	. Tern	nék áttekintés	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.	Csatlakozások	46
2. Sz	állítás	és telepítés	49
2. 1	. Kics	omagolás	49
		A szállítmány tartalma	49
2. 2	. Tele	pítési hely	50
2. 3	. Szá	lítás	52
		A rázógép kezelése	52
2. 4	. Vízs	zintezés	53
2. 5	. Háló	ozati csatlakozás	53
2. 6	. Alap) beállítás	54
2.7	. Táro	lás	59
2. 8	. Elki	ildés	59
2. 8 3. Ke	. Elkü zelés.	ildés	59 60
2. 8 3. Ke 3. 1	zelés. . Be-/	ildés	59 60 60
2. 8 3. Ke 3. 1 3. 2	zelés. zelés. . Be-/ . Grat	ildés	59 60 60 60
2.8 3. Ke 3. 1 3. 2	zelés. zelés. . Be-/ . Grat 3. 2. 1.	ildés Kikapcsolás ïkus felhasználói felület A rázógép legfontosabb üzemi paramétereinek beállítása	59 60 60 60 61
2.8 3. Ke 3. 1 3. 2	zelés. zelés. Be-/ Grat 3. 2. 1. 3. 2. 2.	ildés Kikapcsolás iikus felhasználói felület A rázógép legfontosabb üzemi paramétereinek beállítása Üzemállapot	59 60 60 61 66
2.8 3. Ke 3.1 3.2	 Elki zelés. Be-/ Grat 3. 2. 1. 3. 2. 2. 3. 2. 3. 	ildés Kikapcsolás ikus felhasználói felület A rázógép legfontosabb üzemi paramétereinek beállítása Üzemállapot Beállítások	59 60 60 61 66 68
2.8 3. Ke 3. 1 3.2	 Elki zelés. Be-/ Grat 3. 2. 1. 3. 2. 2. 3. 2. 3. 3. 2. 4. 	ildés Kikapcsolás. Fikus felhasználói felület A rázógép legfontosabb üzemi paramétereinek beállítása Üzemállapot Beállítások Programok	59 60 60 61 66 68 78
2. 8 3. Ke 3. 1 3. 2 3. 3	zelés. Be-/ Grat 3. 2. 1. 3. 2. 2. 3. 2. 3. 3. 2. 4. . Tart	Kikapcsolás . Fikus felhasználói felület A rázógép legfontosabb üzemi paramétereinek beállítása Üzemállapot Beállítások Programok ozékok	59 60 60 61 66 68 78 88
2. 8 3. Ke 3. 1 3. 2 3. 3	 Elki zelés. Be-/ Grat 3. 2. 1. 3. 2. 2. 3. 2. 3. 3. 2. 4. Tart 3. 3. 1. 	Kikapcsolás . Kikapcsolás . Fikus felhasználói felület A rázógép legfontosabb üzemi paramétereinek beállítása Üzemállapot Beállítások Programok Ozékok A tálca felszerelése.	59 60 60 61 66 68 78 88 89
2.8 3. Ke 3.1 3.2 3.3	 Elki zelés. Be-/ Grat 3. 2. 1. 3. 2. 2. 3. 2. 3. 3. 2. 4. Tart 3. 3. 1. 3. 3. 2. 	Kikapcsolás. Fikus felhasználói felület A rázógép legfontosabb üzemi paramétereinek beállítása Üzemállapot Beállítások Programok ozékok A tálca felszerelése. A dugattyútartó és az edény behelyezése	59 60 60 61 66 68 78 88 89 92
2. 8 3. Ke 3. 1 3. 2 3. 3	 Elki zelés. Be-/ Graf 3. 2. 1. 3. 2. 2. 3. 2. 3. 3. 2. 4. Tart 3. 3. 1. 3. 3. 2. 3. 3. 2. 3. 3. 3. 	Kikapcsolás . Kikapcsolás . Kikus felhasználói felület A rázógép legfontosabb üzemi paramétereinek beállítása Üzemállapot Beállítások Programok Ozékok A tálca felszerelése. A dugattyútartó és az edény behelyezése Szögletes lombiktartó felszerelése.	59 60 60 61 66 68 78 88 89 92 92
2.8 3. Ke 3.1 3.2 3.3	 Elki zelés. Be-/ Grat 3. 2. 1. 3. 2. 2. 3. 2. 3. 3. 2. 4. Tart 3. 3. 1. 3. 3. 2. 3. 3. 3. 3. 3. 3. 3. 3. 4. 	Kikapcsolás. řikus felhasználói felület A rázógép legfontosabb üzemi paramétereinek beállítása Üzemállapot Beállítások Programok Ozékok A tálca felszerelése. A dugattyútartó és az edény behelyezése Szögletes lombiktartó felszerelése. A kémcsőállvány felszerelése.	59 60 60 61 66 68 78 88 89 92 94 94 96
2. 8 3. Ke 3. 1 3. 2 3. 3	 Elki zelés. Be-/ Grat 3. 2. 1. 3. 2. 2. 3. 2. 3. 3. 2. 4. Tart 3. 3. 1. 3. 3. 1. 3. 3. 2. 3. 3. 3. 3. 3. 4. 3. 3. 5. 	Kikapcsolás. Fikus felhasználói felület A rázógép legfontosabb üzemi paramétereinek beállítása Üzemállapot Beállítások Programok ozékok A tálca felszerelése. A dugattyútartó és az edény behelyezése Szögletes lombiktartó felszerelése. A kémcsőállvány felszerelése. Állítható szögű kémcsőállvány-tartó felszerelése	59 60 60 61 66 68 78 88 89 92 94 92 94 96 97
2.8 3. Ke 3.1 3.2 3.3	 Elki zelés. Be-/ Grat 3. 2. 1. 3. 2. 2. 3. 2. 3. 3. 2. 4. 3. 3. 1. 3. 3. 1. 3. 3. 2. 3. 3. 3. 3. 3. 4. 3. 3. 5. 3. 3. 6. 	Kikapcsolás. řikus felhasználói felület A rázógép legfontosabb üzemi paramétereinek beállítása Üzemállapot Beállítások Programok Ozékok A tálca felszerelése. A dugattyútartó és az edény behelyezése Szögletes lombiktartó felszerelése. A kémcsőállvány felszerelése. Állítható szögű kémcsőállvány-tartó felszerelése Lombiktartó felszerelése Mikrotiter/Deepwell tálcákhoz	59 60 61 66 68 78 88 89 92 94 96 97 98
2. 8 3. Ke 3. 1 3. 2 3. 3	 Elki zelés. Be-/ Graf 3. 2. 1. 3. 2. 2. 3. 2. 3. 3. 2. 4. Tart 3. 3. 1. 3. 3. 1. 3. 3. 2. 3. 3. 4. 3. 3. 4. 3. 3. 5. 3. 3. 6. 3. 3. 7. 	Kikapcsolás. Fikus felhasználói felület A rázógép legfontosabb üzemi paramétereinek beállítása Üzemállapot Beállítások Programok ozékok A tálca felszerelése. A dugattyútartó és az edény behelyezése Szögletes lombiktartó felszerelése. A kémcsőállvány felszerelése. Állítható szögű kémcsőállvány-tartó felszerelése Lombiktartó felszerelése	59 60 61 66 68 78 88 89 92 94 96 97 98 99

3. 3. 9. A szétválasztó tölcsértartó felszerelése 1	02
3. 3. 10. Többfunkciós tároló felszerelése 1	05
3. 3. 11. Az öntapadós szőnyeg lerakása és használata 1	06
3. 3. 12. Gázelosztó felszerelése 1	13
3. 4. Feltöltés és rendeltetésszerű használat 1	15
Betöltés	16
Rendeltetésszerű használat1	17
4. Karbantartás és ápolás	21
4. 1. Alapok	22
A tartozékok ellenőrzése	23
4. 2. Tisztítás 1	23
Kezelőfelület 1	23
4. 3. Fertőtlenítés 1	24
4. 4. Dekontaminálás 1	24
4. 5. Autoklávozás 1	25
4. 6. Hőmérsékletkalibrálás 1	26
4. 6. 1. Bevált hőmérsékletkalibrálási eljárások	26
4. 6. 2. A hőmérsékletkalibrálás végrehajtása	26
4. 6. 3. A hőmérsékletkalibrálás lekérdezése 1	30
4. 7. A készülékszoftver telepítése 1	30
4. 8. Tálcacsere	34
4. 9. Szerviz	34
4. 10. Elküldés és megsemmisítés	34
5. A NIDAK OKAI ES Elharitasuk 13	35
GPL (General Public License, nyilvános licenc) 13	37
Tárgymutató13	38

Előszó

A kézikönyvről

A kezelési útmutató a következő rázógépeket és tartozékaikat ismerteti (lásd "Tartozékok" a 17. oldalon):

Rázó	Cikkszám
Thermo Scientific Solaris 2000 rázó, 100–240 V ±10 %, 50 / 60 Hz	SK2000
Thermo Scientific Solaris 4000 rázó, 100–240 V ±10 %, 50 / 60 Hz	SK4000
Thermo Scientific Solaris 2000 l rázó, 100–120, 200–240 V ±10 %, 50 / 60 Hz	SK2001
Thermo Scientific Solaris 4000 l rázó, 100–120, 200–240 V ±10 %, 50 / 60 Hz	SK4001
Thermo Scientific Solaris 2000 R rázó, 100–240 V ±10 %, 50 / 60 Hz	SK2002
Thermo Scientific Solaris 4000 R rázó, 100–240 V ±10 %, 50 / 60 Hz	SK4002

Az ebben az útmutatóban található leírások és utasítások nem a teljes terméknévre, hanem az egyes rázógépekre vonatkoznak, osztályonként:

- Nyitott rázógép
- Inkubált rázógép (csak fűtött)
- Hűtött rázógép (hűtött és fűtött)

Rendeltetésszerű használat

Nyitott rázógép

Ezt a készüléket kutatásban és gyártásban használják, oldatok 2D körpályán történő rázására és egyidejűleg a fordulatszám és az idő vezérlésére.

Inkubált rázógép (csak fűtött) és hűtött rázógép (hűtött és fűtött)

Ezt a készüléket a kutatásban és a gyártásban használják oldatok kétdimenziós körpályán történő rázására, és ezzel egyidejűleg a fordulatszám, az időtartam és a hőmérséklet vezérlésére.

Szükséges tudnivalók

A rázót kizárólag megfelelően képzett személyek kezelhetik.

Jelek és kifejezések

Jelző kifejezés	Veszélyfokozat
FIGYELMEZTETÉS	Olyan veszélyhelyzetre utal, ami halált vagy súlyos sérülést okozhat, ha nem kerülik el.
VIGYÁZAT	Olyan veszélyhelyzetre utal, ami könnyű vagy közepes sérülést okozhat, ha nem kerülik el.
MEGJEGYZÉS	Fontos, nem veszélyhez kapcsolódó információkra hívja fel a figyelmet.

A rázón és tartozékain elhelyezett jelek

Tartsa be jelen útmutató előírásait, hogy ne veszélyeztesse saját magát és a környezetét.

8	További információt a kezelési útmutatóban talál
	Húzza ki a hálózati csatlakozót
	Általános veszélyek
	A forró felületek égési sérülést okozhatnak
	Ujj- vagy kézsérülések veszélye a mechanikus alkatrészek véletlenszerű záródása következtében

A kezelési útmutatóban használt jelek

Tartsa be az útmutató előírásait, hogy ne veszélyeztesse a környezetét.

Általános veszélyek	4	Áramütés veszélye
Biológiai veszélyeztetés		Vágási sérülések veszélye
Éghető anyagok veszélye	i	Fontos, nem veszélyhez kapcsolódó információkra hívja fel a figyelmet.
Ujj- vagy kézsérülések veszélye a mechanikus alkatrészek véletlenszerű záródása következtében.		

Biztonsági előírások



Ezen biztonsági előírások figyelmen kívül hagyása súlyos sérüléssel vagy halállal járó veszélyes helyzetet okozhat, ha nem kerülik el.

FIGYELMEZTETÉS

- Tartsa be a biztonsági előírásokat.
- A rázót csak rendeltetésszerűen szabad használni. A rendeltetési céltól eltérő használat károkat, kontaminációt és halálos sérüléseket okozhat.
- A rázót kizárólag megfelelően képzett személyek kezelhetik.
- Az üzemeltető kötelessége biztosítani a megfelelő személyi védőfelszerelés viselését. Vegye figyelembe a WHO "Laboratory Biosafety Manual" és saját országának előírásait.



Hibás áramellátás miatti meghibásodás.

Biztosítsa, hogy a rázót csak megfelelően földelt dugaszoló aljzathoz csatlakoztassák.

FIGYELMEZTETÉS



Veszélyes anyagok kezelése okozta veszély.

Gondosan tisztítsa meg a rázót és a tartozékokat ha korrozív mintákkal (sóoldat, sav, lúg) dolgozik.

FIGYELMEZTETÉS

- A rázó nem inertizált vagy robbanásbiztos. Soha ne használja a rázót robbanásveszélyes környezetben.
- Ne rázzon toxikus vagy radioaktív anyagokat ill. patogén mikroorganizmusokat megfelelő biztonsági intézkedések nélkül.
- Amennyiben veszélyes anyagot ráz, vegye figyelembe a "Laboratory Biosafety manual" kézikönyvet (a WHO kiadványa) és a rá vonatkozó nemzeti ajánlásokat. Ha a WHO "Laboratory Biosafety Manual" szerint II-es kockázati csoportba tartozó mikrobiológiai mintákat ráz, használjon aeroszoltömör biotömítéseket. A "Laboratory Biosafety Manual" a WHO internetes portálján érhető el <u>(www.who.int)</u>. A magas rizikócsoportba tartozó anyagok esetén hozzon további biztonsági intézkedéseket.
- Ha a rázógépet vagy annak részeit toxikus vagy patogén anyagok szennyezték be, megfelelő fertőtlenítési és szennyezésmentesítési intézkedéseket kell végrehajtani (lásd "Dekontaminálás" a 124. oldalon; "Fertőtlenítés" a 124. oldalon).
- Veszélyhelyzet kialakulása esetén szakítsa meg a rázó áramellátását és azonnal hagyja el a területet.



FIGYELMEZTETÉS

Fertőző anyagok okozta egészségkárosodás.

Ha véletlenül folyadékok vagy más anyagok kerülnek a tálca alá, azonnal kapcsolja ki a rázógépet, válassza le a hálózatról és távolítsa el a tálcát (lásd "Tálcák" a 18. oldalon).

Törölje fel a kiömlött anyagokat a bevett laboratóriumi eljárások szerint. Viseljen megfelelő személyi védőfelszerelést.



FIGYELMEZTETÉS

A rázott éghető vagy robbanásveszélyes anyagok vagy közegek egészségkárosodást okozhatnak.

Ne rázzon éghető vagy robbanásveszélyes anyagokat vagy közegeket.



Forró felületek okozta égési sérülések.

VIGYÁZAT

A hőmérséklet-szabályozott rázógépben rendkívül felforrósodhatnak az olyan felületek, mint például a tálca és a burkolat alatti tartozékok. Ne érintse meg a forró felületeket! Várja meg, amíg lehűlnek a forró felületek.



Az üvegcserepek vágott sérüléseket okozhatnak.

VIGYÁZAT

A meglazult tálca és tartozékok, pl. lombiktartó, az üvegek töréséhez vezethetnek, ha azok kiesnek a rázóból.

Biztosítsa, hogy a tálca és a tartozékok megfelelően legyenek felszerelve, a megfelelő szerszámok és csavarok használatával. Biztosítsa, hogy a lombiktartó alkalmas legyen az edény és a minta súlyának felvételére a kívánt fordulatszámon. Az egyedi töltetű lombiktartók esetleg csak olyan fordulatszámra alkalmasak, ami a rázó stabilitási határértéke alatt van. További információkért lásd a következő szakaszt: "Feltöltés és rendeltetésszerű használat" a 115. oldalon.

Figyeljen a szokatlan zajokra. Ezek a tálca és a tartozékok meglazulására utalhatnak.



A rázógép károsodása vagy meghibásodása károsodott kezelőpult következtében.

- VIGYÁZAT
- Ne működtesse a rázógépet, ha károsodott a kezelőpult!
- Kapcsolja ki a rázót. Húzza ki a hálózati csatlakozót. A kezelőpultot arra feljogosított szerviztechnikussal cseréltesse ki.



A törött vagy nem záró edények biológiai veszélyt jelentenek.

A szakszerűtlenül beszerelt tartozékok a minták kiömlését okozhatják.

- Biztosítsa, hogy a tartozékok megfelelően legyenek felszerelve, a megfelelő szerszámok és csavarok használatával.
- Biztosítsa, hogy a tartozékok megfelelően rögzüljenek a tálcára.
- Az edényeket mindig csak az adott méretnek megfelelő tartozékokkal együtt használja.
- Az edények legyenek épek és megfelelően felszerelve.



VIGYÁZAT

A szakszerűtlen feltöltés és a sérült tartozékok befolyásolhatják a biztonságot.

- Mindig ügyeljen arra, hogy a terhelés (tartozékok és minták) a lehető legegyenletesebben legyen elosztva, különösen kétlépcsős tálca üzemeltetése esetén.
 - Ne használjon olyan tartozékot, amin korrózió nyoma vagy repedés látszik. További információkért forduljon ügyfélszolgálatunkhoz.
 - Kizárólag akkor használja a rázógépet, ha azt megfelelően töltötték fel.
 - Soha ne terhelje túl a rázót.
 - Mielőtt a rázót használatba veszi, győződjön meg arról, hogy megfelelően szerelték fel a tartozékokat. A "Tartozékok" a 88. oldalon utasításait be kell tartani.

VIGYÁZAT



Az üzemeltetés alapfeltételeinek figyelmen kívül hagyása sérülést okozhat.

VIGYÁZAT

- Soha ne használja a rázót szabályosan felszerelt tálca nélkül.
- Soha ne használja a rázót, ha a burkolat részei sérültek vagy azokat leszerelték.
- Soha ne használjon meghibásodott burkolatrugóval rendelkező rázógépet! A meghibásodott burkolatrugó nem alkalmas arra, hogy teljesen nyitott helyzetben biztonságosan megtartsa a burkolatot.
- Működés közben ne mozgassa a rázógépet!
- Ne támaszkodjon a rázóra.
- A rázógépet csak akkor töltse fel vagy ürítse ki, ha az teljesen megállt, és a megállást a kezelőpult is megerősíti.
- Működés közben ne tegyen semmit a rázógépre!
- Működés közben ne érintse meg a rázógép tálcáját vagy tartozékait!
- A felhasználó nem nyithatja ki a rázó házát.



A nem kompatibilis tartozékok befolyásolhatják a biztonságot.

A rázóhoz kizárólag a Thermo Fisher Scientific által jóváhagyott tartozékokat használjon. Az aktuális listákat az alábbi címen érheti el: <u>http://thermofisher.com</u>.

MEGJEGYZÉS



A rázó kikapcsolásához:

Nyomja meg a STOP gombot. A rázót a főkapcsolójával kapcsolja ki. Húzza ki a hálózati csatlakozót. Veszély esetén szakítsa meg az áramellátást.

MEGJEGYZÉS

1.1. Műszaki adatok

Thermo Scientific Solaris 2000 rázó					
Fordulatszám tartomány	15-525 ford./perc	l			
Működési idő	99 óra 59 perc (1 perces lép	oésekben) vagy tartós üzem			
Zajszint max. fordulatszámnál	50 dB (A) (1 m a készülék e	lőtt, 1,6 m magasságban)			
Max. terhelés (tálcával, tartozékokkal és mintákkal)	25 kg (55 lbs)				
Elektromos csatlakozás	100-240 V ±10 %, 50 / 60 k	Ηz			
Teljesítményfelvétel	60 W				
Környezeti feltételek					
Tárolás és szállítás esetében	Hőmérséklet: -10 °C - 55 °(Légnedvesség: 15 - 85 %	2			
Használat közben	Beltéri használatra Tengerszint feletti 3000 m magasságig Hőmérséklet: 5 °C - 40 °C Maximális relatív légnedvesség 80 %, 31 °C-ig, lineárisan csökken 50 % relatív légnedvességig, 40 °C-nál.				
Szennyeződési fok	2				
Túlfeszültségi kategória	Ш				
IP	20				
Interfészek	USB	Ethernet			
	2x USB-A 2.0	RJ45			
Méretek	Hossz	Szélesség	Magasság		
	47 cm (18,5 in)	37 cm (14,5 in)	15 cm (5,5 in)		
Súly	20,9 kg (46,0 lbs)				

1. táblázat: Műszaki adatok Solaris 2000

Thermo Scientific Solaris 4000 rázó

Fordulatszám tartomány	15-525 ford./perc		
Működési idő	99 óra 59 perc (1 perces lép	oésekben) vagy tartós üzem	
Zajszint max. fordulatszámnál	54 dB (A) (1 m a készülék e	lőtt, 1,6 m magasságban)	
Max. terhelés (tálcával, tartozékokkal és mintákkal)	43 kg (95 lbs)		
Elektromos csatlakozás	100–240 V ±10 %, 50 / 60 H	Ηz	
Teljesítményfelvétel	80 W		
Környezeti feltételek			
Tárolás és szállítás esetében	Hőmérséklet: -10 °C - 55 °C Légnedvesség: 15 - 85 %	2	
Használat közben	Beltéri használatra Tengerszint feletti 3000 m r Hőmérséklet: 5 °C - 40 °C Maximális relatív légnedves légnedvességig, 40 °C-nál.	nagasságig ség 80 %, 31 °C-ig, lineárisan	ı csökken 50 % relatív
Szennyeződési fok	2		
Túlfeszültségi kategória	Ш		
IP	20		
Interfészek	USB	Ethernet	
	2x USB-A 2.0	RJ45	
Méretek	Hossz	Szélesség	Magasság
	65 cm (25,6 in)	58 cm (22,8 in)	18 cm (7 in)
Súly	75,1 kg (165,5 lbs)		

2. táblázat: Műszaki adatok Solaris 4000

Thermo Scientific Solaris 2000 I rázó

Fordulatszám tartomány	15-525 ford./perc		
Beállítható hőmérsékleti tartomány	30-60 °C		
Hőmérséklet-stabilitás a lombikban 37 °C-nál	Környezeti hőmérséklet 23 °C; 1 órán át stabil hőmérsékletnél a mintakamrában \pm 0,1 °C		
A hőmérséklet egyenletessége a lombikban 37 °C-nál	Környezeti hőmérséklet 23 ° ± 0,5 °C	°C; 1 órán át stabil hőmérsékle	etnél a mintakamrában
Működési idő	99 óra 59 perc (1 perces lép	ésekben) vagy tartós üzem	
Zajszint max. fordulatszámnál	52 dB (A) (1 m a készülék e	lőtt, 1,6 m magasságban)	
Max. terhelés (tálcával, tartozékokkal és mintákkal)	16 kg (35 font), beleértve a 11 x 14 méretű é	és 2,3 kg (5,1 font) tömegű tálo	cát
Elektromos csatlakozás	100–120, 200–240 V ±10 %	, 50 / 60 Hz	
Teljesítményfelvétel	900 W		
Környezeti feltételek			
Tárolás és szállítás esetében	Hőmérséklet: -10 °C - 55 °C Légnedvesség: 15 - 85 %	:	
Használat közben	Beltéri használatra Tengerszint feletti 3000 m magasságig Hőmérséklet: 5 °C - 40 °C Maximális relatív légnedvesség 80 %, 31 °C-ig, lineárisan csökken 50 % relatív légnedvességig, 40 °C-nál.		
Szennyeződési fok	2		
Túlfeszültségi kategória	II		
IP	20		
Interfészek	USB	Ethernet	
	2x USB-A 2.0	RJ45	
Méretek	Hossz	Szélesség	Magasság
	70 cm (27,6 in)	36 cm (14,2 in)	46 cm (18,1 in)
Nyitott burkolattal			79 cm (30,8 in)
Súly	44,9 kg (99 lbs)		

3. táblázat: Műszaki adatok Solaris 2000 l

Thermo Scientific Solaris 4000 I rázó

Fordulatszám tartomány	15-525 ford./perc		
Beállítható hőmérsékleti tartomány	30–60 °C		
Hőmérséklet-stabilitás a lombikban 37 °C-nál	Környezeti hőmérséklet 23 °C; 1 órán át stabil hőmérsékletnél a mintakamrában \pm 0,1 °C		
A hőmérséklet egyenletessége a lombikban 37 °C-nál	Környezeti hőmérséklet 23 ° ± 0,5 °C	C; 1 órán át stabil hőmérsékle	tnél a mintakamrában
Működési idő	99 óra 59 perc (1 perces lép	ésekben) vagy tartós üzem	
Zajszint max. fordulatszámnál	52 dB (A) (1 m a készülék el	őtt, 1,6 m magasságban)	
Max. terhelés (tálcával, tartozékokkal és mintákkal)	32 kg (70,5 font), beleértve a 18 x 18 méretű é	às 4,3 kg (9,5 font) tömegű tálo	cát
Elektromos csatlakozás	100-120, 200-240 V ±10 %	, 50 / 60 Hz	
Teljesítményfelvétel	900 W		
Környezeti feltételek			
Tárolás és szállítás esetében	Hőmérséklet: -10 °C - 55 °C Légnedvesség: 15 - 85 %		
Használat közben	Beltéri használatra Tengerszint feletti 3000 m magasságig Hőmérséklet: 5 °C - 40 °C Maximális relatív légnedvesség 80 %, 31 °C-ig, lineárisan csökken 50 % relatív légnedvességig, 40 °C-nál.		
Szennyeződési fok	2		
Túlfeszültségi kategória	II		
IP	20		
Interfészek	USB	Ethernet	
	2x USB-A 2.0	RJ45	
Méretek	Hossz	Szélesség	Magasság
	77 cm (30,3 in)	57 cm (22,4 in)	55 cm (21,7 in)
Nyitott burkolattal			94 cm (36,7 in)
Súly	69.1 kg (150.1 lbg)		

4. táblázat: Műszaki adatok Solaris 4000 l

Thermo Scientific Solaris 2000 R rázó

Fordulatszám tartomány	15-525 ford./perc		
Beállítható hőmérsékleti tartomány	5–60 °C		
Hőmérséklet-stabilitás a lombikban 37 °C-nál	Környezeti hőmérséklet 23 °C; 1 órán át stabil hőmérsékletnél a mintakamrában \pm 0,1 °C		
A hőmérséklet egyenletessége a lombikban 37 °C-nál	Környezeti hőmérséklet 23 ° ± 0,5 °C	°C; 1 órán át stabil hőmérsékle	tnél a mintakamrában
Működési idő	99 óra 59 perc (1 perces lép	ésekben) vagy tartós üzem	
Zajszint max. fordulatszámnál	56 dB (A) (1 m a készülék e	lőtt, 1,6 m magasságban)	
Max. terhelés (tálcával, tartozékokkal és mintákkal)	16 kg (35 font), beleértve a 11 x 14 méretű é	és 2,3 kg (5,1 font) tömegű tálo	sát
Elektromos csatlakozás	100-240 V ±10 %, 50 / 60 H	łz	
Teljesítményfelvétel	350 W		
Környezeti feltételek			
Tárolás és szállítás esetében	Hőmérséklet: -10 °C - 55 °C Légnedvesség: 15 - 85 %	:	
Használat közben	Beltéri használatra Tengerszint feletti 3000 m magasságig Hőmérséklet: 5 °C - 40 °C Maximális relatív légnedvesség 80 %, 31 °C-ig, lineárisan csökken 50 % relatív légnedvességig, 40 °C-nál.		
Szennyeződési fok	2		
Túlfeszültségi kategória	II		
IP	20		
Interfészek	USB	Ethernet	
	2x USB-A 2.0	RJ45	
Méretek	Hossz	Szélesség	Magasság
	70 cm (27,6 in)	36 cm (14,2 in)	46 cm (18,1 in)
Nyitott burkolattal			79 cm (30,8 in)
Nyitott burkolattal Súly	47,2 kg (104,1 lbs)		79 cm (30,8 in)

5. táblázat: Műszaki adatok Solaris 2000 R

Thermo Scientific Solaris 4000 R rázó

Fordulatszám tartomány	15-525 ford./perc			
Beállítható hőmérsékleti tartomány	4–60 °C			
Hőmérséklet-stabilitás a lombikban 37 °C-nál	Környezeti hőmérséklet 23 °(± 0,1 °C	C; 1 órán át stabil hőmérséklel	nél a mintakamrában	
A hőmérséklet egyenletessége a lombikban 37 °C-nál	Környezeti hőmérséklet 23 °C; 1 órán át stabil hőmérsékletnél a mintakamrában \pm 0,5 °C			
Működési idő	99 óra 59 perc (1 perces lépésekben) vagy tartós üzem			
Zajszint max. fordulatszámnál	56 dB (A) (1 m a készülék elő	itt, 1,6 m magasságban)		
Max. terhelés (tálcával, tartozékokkal és mintákkal)	32 kg (70,5 font), beleértve a 18 x 18 méretű és 4,3 kg (9,5 font) tömegű tálcát			
Elektromos csatlakozás	100–240 V ±10 %, 50 / 60 Hz			
Teljesítményfelvétel	750 W			
Környezeti feltételek				
Tárolás és szállítás esetében	Hőmérséklet: -10 °C - 55 °C Légnedvesség: 15 - 85 %			
Használat közben	Beltéri használatra Tengerszint feletti 3000 m magasságig Hőmérséklet: 5 °C - 40 °C Maximális relatív légnedvesség 80 %, 31 °C-ig, lineárisan csökken 50 % relatív légnedvességig, 40 °C-nál.			
Szennyeződési fok	2			
Túlfeszültségi kategória	II			
IP	20			
Interfészek	USB	Ethernet		
	2x USB-A 2.0	RJ45		
Méretek	Hossz	Szélesség	Magasság	
	77 cm (30,3 in)	57 cm (22,4 in)	55 cm (21,7 in)	
Nyitott burkolattal			94 cm (36,7 in)	
Súly	74,9 kg (165,1 lbs)			

6. táblázat: Műszaki adatok Solaris 4000 R

1.2. Tartozékok



A nem kompatibilis tartozékok befolyásolhatják a biztonságot. A rázóhoz kizárólag a Thermo Fisher Scientific által jóváhagyott tartozékokat használjon.

VIGYÁZAT

Az aktuális listákat az alábbi címen érheti el: http://thermofisher.com.

1. 2. 1. Tálcák

Tálca		
Solaris 2000	Súly	
Univerzális tálca, Thermo Scientific Solaris 12x14	2,7 kg (5,9 lbs)	SK1214
Kétlépcsős univerzális tálca, Thermo Scientific Solaris 12x14	6,4 kg (14,1 lbs)	SK1214D
Kiegészítő készlet kétlépcsős univerzális tálcához, Thermo Scientific Solaris 12x14		SK1214DK
Univerzális tálca, Thermo Scientific Solaris 18x18	4,3 kg (9,4 lbs)	SK1818
Kétlépcsős univerzális tálca, Thermo Scientific Solaris 18x18	9,7 kg (21,3 lbs)	SK1818D
Kiegészítő készlet kétlépcsős univerzális tálcához, Thermo Scientific Solaris 18x18		SK1818DK
Univerzális tálca, Thermo Scientific Solaris 18x24	6,0 kg (13,2 lbs)	SK1824
Solaris 4000		
Univerzális tálca, Thermo Scientific Solaris 18x30	7,5 kg (16,5 lbs)	SK1830
Kétlépcsős univerzális tálca, Thermo Scientific Solaris 18x30	16,3 kg (35,9 lbs)	SK1830D
Kiegészítő készlet kétlépcsős univerzális tálcához, Thermo Scientific Solaris 18x30		SK1830DK
Univerzális tálca, Thermo Scientific Solaris 36x24	12,0 kg (26,4 lbs)	SK3624
Solaris 2000 I / 2000 R		
Univerzális tálca, Thermo Scientific Solaris 11x14	2,3 kg (5,1 lbs)	SK1114
Solaris 4000 1 / 4000 R		
Univerzális tálca, Thermo Scientific Solaris 18x18	4,3 kg (9,5 lbs)	SK1818
Utólag felszerelhető készletek és tartozékok		
Lombiktartó pótalkatrész készlet (csavarok)		SK1001
Pótalkatrészkészlet tálcához, kicsi (tálcacsavarok, szerszám)		SK0100
Pótalkatrészkészlet tálcához, nagy (tálcacsavarok, szerszám)		SK0101
Csavarhúzó a tartozékokhoz		75004131

7. táblázat: Az elérhető tálcák

1. 2. 2. Lombiktartó



Solaris 2000 tálcákhoz

	Cikkszám	Univerzális			Kétlép	csős
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
10 ml Erlenmeyer lombik(ok)	30150BI	72	113	157	140	226
25 ml Erlenmeyer lombik(ok)	30151	42	64	80	80	124
50 ml Erlenmeyer lombik(ok)	30152BI	42	64	80	80	124
125 ml Erlenmeyer lombik(ok)	30153	15	32	40	30	52
250 ml Erlenmeyer lombik(ok)	30154BI	9	16	24	16	32
300 ml Erlenmeyer lombik(ok)	30155	9	16	20	16	32
500 ml Erlenmeyer lombik(ok)	30156BI	9	16	20	16	32
1 I Erlenmeyer lombik(ok)	30157BI	4	9	10	8	16
2 I Erlenmeyer lombik(ok)	30158	3	5	6	-	-
4 I Erlenmeyer lombik(ok)	30159	1	4	4	-	-
5 I Erlenmeyer lombik(ok)	30159B	1	2	4	-	-
6 I Erlenmeyer lombik(ok)	30160	1	2	2	-	-
2800 ml Fernbach lombik(ok)	30162	1	4	4	-	-
2500 ml tenyésztő lombik(ok), alacsony kivitel	30161	1	1	2	-	-

8. táblázat: A Solaris 2000 tálcákhoz elérhető lombiktartók

Solaris 4000 tálcákhoz

	Cikkszám	Univerzális		Kétlépcsős
		18 x 30	36 x 24	18 x 30
10 ml Erlenmeyer lombik(ok)	30150BI	203	187	402
25 ml Erlenmeyer lombik(ok)	30151	112	187	220
50 ml Erlenmeyer lombik(ok)	30152BI	112	187	220
125 ml Erlenmeyer lombik(ok)	30153	46	83	92
250 ml Erlenmeyer lombik(ok)	30154BI	28	40	56
300 ml Erlenmeyer lombik(ok)	30155	28	40	56
500 ml Erlenmeyer lombik(ok)	30156BI	28	40	56
1 Erlenmeyer lombik(ok)	30157BI	14	20	28
2 Erlenmeyer lombik(ok)	30158	6	11	12
4 I Erlenmeyer lombik(ok)	30159	6	8	-
5 I Erlenmeyer lombik(ok)	30159B	5	8	-
6 I Erlenmeyer lombik(ok)	30160	3	6	-
2800 ml Fernbach lombik(ok)	30162	6	8	12
2500 ml tenyésztő lombik(ok), alacsony kivitel	30161	3	6	6

9. táblázat: A Solaris 4000 tálcákhoz elérhető lombiktartók

Solaris 2000 I / 2000 R tálcákhoz

	Cikkszám	Univerzális
		11 x 14
10 ml Erlenmeyer lombik(ok)	30150BI	59
25 ml Erlenmeyer lombik(ok)	30151	35
50 ml Erlenmeyer lombik(ok)	30152BI	35
125 ml Erlenmeyer lombik(ok)	30153	15
250 ml Erlenmeyer lombik(ok)	30154BI	7
300 ml Erlenmeyer lombik(ok)	30155	7
500 ml Erlenmeyer lombik(ok)	30156BI	7
1 Erlenmeyer lombik(ok)	30157BI	4
2 I Erlenmeyer lombik(ok)	30158	2
4 I Erlenmeyer lombik(ok)	30159	-
5 I Erlenmeyer lombik(ok)	30159B	-
6 I Erlenmeyer lombik(ok)	30160	-
2800 ml Fernbach lombik(ok)	30162	1
2500 ml tenyésztő lombik(ok), alacsony kivitel	30161	1

10. táblázat: A Solaris 2000 I / 2000 R tálcákhoz elérhető lombiktartók

Solaris 4000 I / 4000 R tálcákhoz

	Cikkszám	Univerzális
		18 x 18
10 ml Erlenmeyer lombik(ok)	30150BI	113
25 ml Erlenmeyer lombik(ok)	30151	64
50 ml Erlenmeyer lombik(ok)	30152BI	64
125 ml Erlenmeyer lombik(ok)	30153	32
250 ml Erlenmeyer lombik(ok)	30154BI	16
300 ml Erlenmeyer lombik(ok)	30155	16
500 ml Erlenmeyer lombik(ok)	30156BI	16
1 I Erlenmeyer lombik(ok)	30157BI	9
2 I Erlenmeyer lombik(ok)	30158	6
4 I Erlenmeyer lombik(ok)	30159	-
5 I Erlenmeyer lombik(ok)	30159B	-
6 I Erlenmeyer lombik(ok)	30160	-
2800 ml Fernbach lombik(ok)	30162	4
2500 ml tenyésztő lombik(ok), alacsony kivitel	30161	1

11. táblázat: A Solaris 4000 I / 4000 R tálcákhoz elérhető lombiktartók

1. 2. 3. Lombiktartó Mikrotiter/Deepwell tálcákhoz



Solaris 2000 tálcákhoz

	Cikkszám	Univerzális		Kétlépcsős		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Mikrotiter- / Deepwell-tálca	30175	7	12	16	14	24

12. táblázat: A lombiktartók kapacitása Mikrotiter/Deepwell tálcákhoz Solaris 2000 tálcák esetén

Solaris 2000 I / 2000 R tálcákhoz

	Cikkszám	Univerzális
		11 x 14
Mikrotiter- / Deepwell-tálca	30175	6

13. táblázat: Mikrotiter/Deepwell tálcatartók kapacitása Solaris 2000 I/2000 R tálcák esetén

Solaris 4000 tálcákhoz

	Cikkszám	Unive	Univerzális	
		18 x 30	36 x 24	18 x 30
Mikrotiter- / Deepwell-tálca	30175	22	36	42

14. táblázat: Mikrotiter/Deepwell tálcatartók kapacitása Solaris 4000 tálcák esetén

Solaris 4000 I / 4000 R tálcákhoz

	Cikkszám	Univerzális
		18 x 18
Mikrotiter- / Deepwell-tálca	30175	10

15. táblázat: Mikrotiter/Deepwell tálcatartók kapacitása Solaris 4000 I/4000 R tálcák esetén

1. 2. 4. Kémcsőállványok

Solaris 2000 tálcákhoz



	Cikkszám	Univerzális			Kétlépcsős	
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
feles méret						
10–13 mm, piros, 6 x 6-os elrendezés	30181	8	12	15	14	24
14–16 mm, narancs, 6 x 6-os elrendezés	30183	5	9	11	9	17
17–20 mm, fehér, 4 x 5-ös elrendezés	30185	7	11	14	13	21
21–25 mm, kék, 4 x 4-os elrendezés	30187	6	9	11	10	17
26–30 mm, zöld, 3 x 3-os elrendezés	30189	6	9	12	11	18
Mikrotiter kémcsövekhez, 1,5 ml, kék, 4 x 6-os elrendezés	30191	6	10	13	11	19
Méret						
10-13 mm, 6 x 12-es elrendezés	30180BI	3	7	10	6	13
14-16 mm, 6 x 12-es elrendezés	30182	3	4	6	6	8
17-20 mm, 4 x 10-es elrendezés	30184	3	5	7	6	8
21-25 mm, 4 x 10-es elrendezés	30186	2	3	5	4	6
26-30 mm, 3 x 8-es elrendezés	30188	3	4	6	5	8
Mikrotiter kémcsövekhez, 1,5 ml, 8 x 12-es elrendezés	30190	3	5	7	6	9

16. táblázat: A Solaris 2000 tálcákhoz elérhető kémcsőállványok

Solaris 4000 tálcákhoz

	Cikkszám	Univer	Univerzális	
		18 x 30	36 x 24	18 x 30
feles méret				
10–13 mm, piros, 6 x 6-os elrendezés	30181	21	32	40
14–16 mm, narancs, 6 x 6-os elrendezés	30183	14	20	26
17–20 mm, fehér, 4 x 5-ös elrendezés	30185	18	20	34
21–25 mm, kék, 4 x 4-os elrendezés	30187	14	22	25
26–30 mm, zöld, 3 x 3-os elrendezés	30189	15	24	30
Mikrotiter kémcsövekhez, 1,5 ml, kék, 4 x 6-os elrendezés	30191	18	24	34
Méret				
10-13 mm, 6 x 12-es elrendezés	30180BI	14	20	26
14-16 mm, 6 x 12-es elrendezés	30182	9	12	16
17-20 mm, 4 x 10-es elrendezés	30184	9	15	18
21-25 mm, 4 x 10-es elrendezés	30186	7	9	13
26-30 mm, 3 x 8-es elrendezés	30188	7	10	13
Mikrotiter kémcsövekhez, 1,5 ml, 8 x 12-es elrendezés	30190	9	12	17

17. táblázat: A Solaris 4000 tálcákhoz elérhető kémcsőállványok

Solaris 2000 I / 2000 R tálcákhoz

	Cikkszám	Univerzális
		11 x 14
feles méret		
10–13 mm, piros, 6 x 6-os elrendezés	30181	6
14–16 mm, narancs, 6 x 6-os elrendezés	30183	2
17–20 mm, fehér, 4 x 5-ös elrendezés	30185	4
21–25 mm, kék, 4 x 4-os elrendezés	30187	2
26–30 mm, zöld, 3 x 3-os elrendezés	30189	4
Mikrotiter kémcsövekhez, 1,5 ml, kék, 4 x 6-os elrendezés	30191	3
Méret		
10-13 mm, 6 x 12-es elrendezés	30180BI	3
14-16 mm, 6 x 12-es elrendezés	30182	2
17-20 mm, 4 x 10-es elrendezés	30184	2
21-25 mm, 4 x 10-es elrendezés	30186	1
26-30 mm, 3 x 8-es elrendezés	30188	2
Mikrotiter kémcsövekhez, 1,5 ml, 8 x 12-es elrendezés	30190	2

18. táblázat: Solaris 2000 I/2000 R tálcákhoz rendelkezésre álló kémcsőállványok

Solaris 4000 I / 4000 R tálcákhoz

	Cikkszám	Univerzális
		18 x 18
feles méret		
10–13 mm, piros, 6 x 6-os elrendezés	30181	12
14–16 mm, narancs, 6 x 6-os elrendezés	30183	8
17–20 mm, fehér, 4 x 5-ös elrendezés	30185	10
21–25 mm, kék, 4 x 4-os elrendezés	30187	8
26–30 mm, zöld, 3 x 3-os elrendezés	30189	9
Mikrotiter kémcsövekhez, 1,5 ml, kék, 4 x 6-os elrendezés	30191	8
Méret		
10-13 mm, 6 x 12-es elrendezés	30180BI	7
14-16 mm, 6 x 12-es elrendezés	30182	4
17-20 mm, 4 x 10-es elrendezés	30184	5
21-25 mm, 4 x 10-es elrendezés	30186	3
26-30 mm, 3 x 8-es elrendezés	30188	4
Mikrotiter kémcsövekhez, 1,5 ml, 8 x 12-es elrendezés	30190	5

19. táblázat: Solaris 4000 I/4000 R tálcákhoz rendelkezésre álló kémcsőállványok

1. 2. 5. Állítható szögű kémcsőállványok



Solaris 2000 tálcákhoz

	Cikkszám	Univerzális			Kétlépcsős		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18	
10-13 mm, 72 kémcsőhöz	236090	3	4	6	6	8	
16-20 mm, 40 kémcsőhöz	236091	3	4	6	5	8	
21-25 mm, 40 kémcsőhöz	236092	3	3	4	5	6	
26-30 mm, 24 kémcsőhöz	236093	3	3	4	5	6	

20. táblázat: Solaris 2000 tálcákhoz választható állítható szögű kémcsőállványok

Solaris 4000 tálcákhoz

	Cikkszám	Univerzális		Kétlépcsős	
		18 x 30	36 x 24	18 x 30	
10-13 mm, 72 kémcsőhöz	236090	7	14	14	
16-20 mm, 40 kémcsőhöz	236091	7	14	14	
21-25 mm, 40 kémcsőhöz	236092	5	9	10	
26-30 mm, 24 kémcsőhöz	236093	6	9	12	

21. táblázat: Solaris 4000 tálcákhoz választható állítható szögű kémcsőállványok

Solaris 2000 I / 2000 R tálcákhoz

	Cikkszám	Univerzális
		11 x 14
10-13 mm, 72 kémcsőhöz	236090	2
16-20 mm, 40 kémcsőhöz	236091	2
21-25 mm, 40 kémcsőhöz	236092	1
26-30 mm, 24 kémcsőhöz	236093	1

22. táblázat: Solaris 2000 I/2000 R tálcákhoz választható állítható szögű kémcsőállványok

Solaris 4000 I / 4000 R tálcákhoz

	Cikkszám	Univerzális
		18 x 18
10-13 mm, 72 kémcsőhöz	236090	4
16-20 mm, 40 kémcsőhöz	236091	3
21-25 mm, 40 kémcsőhöz	236092	3
26-30 mm, 24 kémcsőhöz	236093	3

23. táblázat: Solaris 4000 I/4000 R tálcákhoz választható állítható szögű kémcsőállványok

1. 2. 6. Állítható edénytartó

Solaris 2000 tálcákhoz

	Cikkszám	Univerzális		Kétlépcsős		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Egy fokozatban állítható magasságú edénytartó 11"x14"	75004104	-	-	-	-	-
Két fokozatban állítható magasságú edénytartó 12"x14"	75004102	1	1	1	-	2
Két fokozatban állítható magasságú edénytartó 18"x18"	75004103	-	1	1	-	-
Állítható egyes edénytartó	75004101	4	4	6	6	8

24. táblázat: Solaris 2000 tálcákhoz választható állítható edénytartók

Solaris 4000 tálcákhoz

	Cikkszám	Univerzális		Kétlépcsős
		18 x 30	36 x 24	18 x 30
Egy fokozatban állítható magasságú edénytartó 11"x14"	75004104	-	-	-
Két fokozatban állítható magasságú edénytartó 12"x14"	75004102	2	2	4
Két fokozatban állítható magasságú edénytartó 18"x18"	75004103	1	2	2
Állítható egyes edénytartó	75004101	8	12	16

25. táblázat: Állítható edénytartók Solaris 4000 tálcákhoz

Solaris 2000 I / 2000 R tálcákhoz

	Cikkszám	Univerzális
		11 x 14
Egy fokozatban állítható magasságú edénytartó 11"x14"	75004104	1
Két fokozatban állítható magasságú edénytartó 12"x14"	75004102	-
Két fokozatban állítható magasságú edénytartó 18"x18"	75004103	-
Állítható egyes edénytartó	75004101	2

26. táblázat: Állítható edénytartók Solaris 2000 I/2000 R tálcákhoz

Solaris 4000 I / 4000 R tálcákhoz

	Cikkszám	Univerzális
		18 x 18
Egy fokozatban állítható magasságú edénytartó 11"x14"	75004104	-
Két fokozatban állítható magasságú edénytartó 12"x14"	75004102	1
Két fokozatban állítható magasságú edénytartó 18"x18"	75004103	1
Állítható egyes edénytartó	75004101	4

27. táblázat: Állítható edénytartók Solaris 4000 I/4000 R tálcákhoz

1. 2. 7. Szögletes edénytartó



Solaris 2000 tálcákhoz

	Cikkszám	Univerzális			Kétlépcsős		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18	
Szögletes palacktartó, 125 ml	75004106	12	16	24	21	32	
Szögletes palacktartó, 250 ml	75004107	9	16	22	18	32	
Szögletes palacktartó, 500 ml	75004108	6	9	12	11	18	
Szögletes palacktartó, 1000 ml	75004109	4	9	11	8	18	

28. táblázat: Solaris 2000 tálcákhoz választható szögletes edénytartók

Solaris 4000 tálcákhoz

	Cikkszám	Unive	rzális	Kétlépcsős
		18 x 30	36 x 24	18 x 30
Szögletes palacktartó, 125 ml	75004106	28	54	56
Szögletes palacktartó, 250 ml	75004107	28	52	56
Szögletes palacktartó, 500 ml	75004108	15	24	30
Szögletes palacktartó, 1000 ml	75004109	15	24	28

29. táblázat: Solaris 4000 tálcákhoz választható szögletes edénytartók

Solaris 2000 I / 2000 R tálcákhoz

	Cikkszám	Univerzális
		11 x 14
Szögletes palacktartó, 125 ml	75004106	9
Szögletes palacktartó, 250 ml	75004107	7
Szögletes palacktartó, 500 ml	75004108	3
Szögletes palacktartó, 1000 ml	75004109	2

30. táblázat: Solaris 2000 I/2000 R tálcákhoz választható szögletes edénytartók

Solaris 4000 I / 4000 R tálcákhoz

	Cikkszám	Univerzális
		18 x 18
Szögletes palacktartó, 125 ml	75004106	16
Szögletes palacktartó, 250 ml	75004107	16
Szögletes palacktartó, 500 ml	75004108	9
Szögletes palacktartó, 1000 ml	75004109	9

31. táblázat: Solaris 4000 I/4000 R tálcákhoz választható szögletes edénytartók

1. 2. 8. Nalgene főzőpohár állványok

Solaris 2000 tálcákhoz



	Cikkszám		Univerzális		Kétlép	ocsős
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Méret 11 x 14/12 x 14			ÁI	lványok/helye	łk	
Nalgene 30 ml-es főzőpohár, 36 hely	75004116	1 / 36	-	-	2/72	-
Nalgene 50 ml-es főzőpohár, 28 hely	75004129	1 / 28	-	-	2 / 56	-
Nalgene 100 ml-es főzőpohár, 16 hely	75004118	1 / 16	-	-	2/32	-
Nalgene 250 ml-es főzőpohár, 9 hely	75004119	1/9	-	-	2/18	-
Nalgene 400 ml-es főzőpohár, 8 hely	75004120	1/8	-	-	2/16	-
Nalgene 600 ml-es főzőpohár, 5 hely	75004121	1/5	-	-	2/10	-
Méret 9 x 18			ÁI	lványok/helye	łk	
Nalgene 30 ml-es főzőpohár, 32 hely	75004110	-	2/64	2 / 64	-	4 / 128
Nalgene 50 ml-es főzőpohár, 24 hely	75004128	-	2 / 48	2 / 48	-	4 / 96
Nalgene 100 ml-es főzőpohár, 15 hely	75004112	-	2/30	2 / 30	-	4 / 60
Nalgene 250 ml-es főzőpohár, 8 hely	75004113	-	2 / 16	2 / 16	-	4 / 32
Nalgene 400 ml-es főzőpohár, 6 hely	75004114	-	2 / 12	2 / 12	-	4 / 24
Nalgene 600 ml-es főzőpohár, 4 hely	75004115	-	2/8	2/8	-	4 / 16

32. táblázat: Solaris 2000 tálcákhoz választható Nalgene főzőpohár állványok

Solaris 4000 tálcákhoz

	Cikkszám	Univerzális		Kétlépcsős	
		18 x 30	36 x 24	18 x 30	
Méret 11 x 14/12 x 14			Állványok/helyek		
Nalgene 30 ml-es főzőpohár, 36 hely	75004116	-	-	-	
Nalgene 50 ml-es főzőpohár, 28 hely	75004129	-	-	-	
Nalgene 100 ml-es főzőpohár, 15 hely	75004118	-	-	-	
Nalgene 250 ml-es főzőpohár, 9 hely	75004119	-	-	-	
Nalgene 400 ml-es főzőpohár, 8 hely	75004120	-	-	-	
Nalgene 600 ml-es főzőpohár, 5 hely	75004121	-	-	-	
Méret 9 x 18			Állványok/helyek		
Nalgene 30 ml-es főzőpohár, 32 hely	75004110	3 / 96	4 / 128	6 / 192	
Nalgene 50 ml-es főzőpohár, 24 hely	75004128	3/72	4 / 96	6 / 144	
Nalgene 100 ml-es főzőpohár, 15 hely	75004112	3/45	4 / 60	6 / 90	
Nalgene 250 ml-es főzőpohár, 8 hely	75004113	3 / 24	4 / 32	6 / 48	
Nalgene 400 ml-es főzőpohár, 6 hely	75004114	3 / 18	4 / 24	6 / 36	
Nalgene 600 ml-es főzőpohár, 4 hely	75004115	3 / 12	4 / 16	6 / 24	

33. táblázat: Solaris 4000 tálcákhoz választható Nalgene főzőpohár állványok

Solaris 2000 I / 2000 R tálcákhoz

	Cikkszám	Univerzális
		11 x 14
Méret 11 x 14/12 x 14		Állványok/helyek
Nalgene 30 ml-es főzőpohár, 36 hely	75004116	1/36
Nalgene 50 ml-es főzőpohár, 28 hely	75004129	1 / 28
Nalgene 100 ml-es főzőpohár, 16 hely	75004118	1 / 16
Nalgene 250 ml-es főzőpohár, 9 hely	75004119	1/9
Nalgene 400 ml-es főzőpohár, 8 hely	75004120	1/8
Nalgene 600 ml-es főzőpohár, 5 hely	75004121	1/5
Méret 9 x 18		Állványok/helyek
Nalgene 30 ml-es főzőpohár, 32 hely	75004110	-
Nalgene 50 ml-es főzőpohár, 24 hely	75004128	-
Nalgene 100 ml-es főzőpohár, 15 hely	75004112	-
Nalgene 250 ml-es főzőpohár, 8 hely	75004113	-
Nalgene 400 ml-es főzőpohár, 6 hely	75004114	-
Nalgene 600 ml-es főzőpohár, 4 hely	75004115	-

34. táblázat: Solaris 2000 I/2000 R tálcákhoz választható Nalgene főzőpohár állványok
Tartozékok

Solaris 4000 I / 4000 R tálcákhoz

	Cikkszám	Univerzális
		18 x 18
Méret 11 x 14/12 x 14		Állványok/helyek
Nalgene 30 ml-es főzőpohár, 36 hely	75004116	-
Nalgene 50 ml-es főzőpohár, 28 hely	75004129	-
Nalgene 100 ml-es főzőpohár, 16 hely	75004118	-
Nalgene 250 ml-es főzőpohár, 9 hely	75004119	-
Nalgene 400 ml-es főzőpohár, 8 hely	75004120	-
Nalgene 600 ml-es főzőpohár, 5 hely	75004121	-
Méret 9 x 18		Állványok/helyek
Nalgene 30 ml-es főzőpohár, 32 hely	75004110	2/64
Nalgene 50 ml-es főzőpohár, 24 hely	75004128	2 / 48
Nalgene 100 ml-es főzőpohár, 15 hely	75004112	2 / 30
Nalgene 250 ml-es főzőpohár, 8 hely	75004113	2 / 16
Nalgene 400 ml-es főzőpohár, 6 hely	75004114	2 / 12
Nalgene 600 ml-es főzőpohár, 4 hely	75004115	2/8

35. táblázat: Solaris 4000 I/4000 R tálcákhoz választható Nalgene főzőpohár állványok

Tartozékok

1. 2. 9. Szétválasztó tölcsértartó

Solaris 2

olaris 2000 tálcákhoz					27	6
	Cikkszám		Kétlépcsős			
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Egyes lombiktartó szétválasztó tölcsérhez, 250 ml és 2000 ml között	75004125	2	3	4	2	5
Szétválasztó tölcsértartó, függőlegesen szerelt		4	4	6	-	-

. Je

36. táblázat: Solaris 2000 tálcákhoz választható szétválasztó tölcsértartók

Solaris 4000 tálcákhoz

	Cikkszám	Univerzális		Kétlépcsős
		18 x 30	36 x 24	18 x 30
Egyes lombiktartó szétválasztó tölcsérhez, 250 ml és 2000 ml között	75004125	5	7	9
Szétválasztó tölcsértartó, függőlegesen szerelt		8	12	-

37. táblázat: Solaris 4000 tálcákhoz választható szétválasztó tölcsértartók

Solaris 2000 I / 2000 R tálcákhoz

	Cikkszám	Univerzális
		11 x 14
Egyes lombiktartó szétválasztó tölcsérhez, 250 ml és 2000 ml között	75004125	-
Szétválasztó tölcsértartó, függőlegesen szerelt		-

38. táblázat: Solaris 2000 I/2000 R tálcákhoz választható szétválasztó tölcsértartók

Solaris 4000 I / 4000 R tálcákhoz

	Cikkszám	Univerzális
		18 x 18
Egyes lombiktartó szétválasztó tölcsérhez, 250 ml és 2000 ml között	75004125	-
Szétválasztó tölcsértartó, függőlegesen szerelt		-

39. táblázat: Solaris 4000 I/4000 R tálcákhoz választható szétválasztó tölcsértartók

1. 2. 10. Többfunkciós tároló



Solaris 2000 tálcákhoz

	Cikkszám	Univerzális			Kétlépcsős		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18	
Többfunkciós tároló 11 x 14/12 x 14	75004123	1	-	-	2	-	
Többfunkciós tároló 9 x 18	75004122	-	2	3	-	4	

40. táblázat: Solaris 2000 tálcákhoz választható többfunkciós tárolók

Solaris 4000 tálcákhoz

	Cikkszám	Univer	zális	Kétlépcsős
		18 x 30	36 x 24	18 x 30
Többfunkciós tároló 11 x 14/12 x 14	75004123	-	-	-
Többfunkciós tároló 9 x 18	75004122	3	6	6

41. táblázat: Solaris 4000 tálcákhoz választható többfunkciós tárolók

Solaris 2000 I / 2000 R tálcákhoz

	Cikkszám	Univerzális
		11 x 14
Többfunkciós tároló 11 x 14/12 x 14	75004123	1
Többfunkciós tároló 9 x 18	75004122	-

42. táblázat: Solaris 2000 I/2000 R tálcákhoz választható többfunkciós tárolók

Solaris 4000 I / 4000 R tálcákhoz

	Cikkszám	Univerzális
		18 x 18
Többfunkciós tároló 11 x 14/12 x 14	75004123	-
Többfunkciós tároló 9 x 18	75004122	2

43. táblázat: Solaris 4000 I/4000 R tálcákhoz választható többfunkciós tárolók

Többfunkciós tárolók tartozékai

	Cikkszám	Univerzális
Tartalék O-gyűrűk (többfunkciós tároló)	75004132	-

44. táblázat: Többfunkciós tárolókhoz választható tartozékok

1. 2. 11. Öntapadós szőnyegek

Általános információk az öntapadós szőnyegek kiválasztásával kapcsolatban:

- A 170 x 280 mm méretű szőnyegek a 11 x 14 és 12 x 14 méretű tálcákhoz ideálisak.
- A 200 x 200 mm méretű szőnyegek a nagyobb, 18 x 18, 18 x 24, 30 x 18 és 36 x 24 méretű tálcákhoz ideálisak.
- Minden szőnyeg tetszőleges méretre vágható.

Solaris 2000 tálcákhoz

	Cikkszám Univerzális K L			Kétlép unive	ocsős, rzális	
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Nagy tapadású szőnyeg, 200 x 200 mm, csomag. egység: 2 db.	75004126	1	4	6	2	8
Nagy tapadású szőnyeg, 280 x 170 mm, csomag. egység: 2 db.	75004127	2	2	4	4	4
Kis tapadású szőnyeg, 200 x 200 mm, csomag. egység: 2 db.	75004111	1	4	6	2	8
Kis tapadású szőnyeg, 280 x 170 mm, csomag. egység: 2 db.	75004117	2	2	4	4	4

A szőnyegek 2 db-os kiszerelésben kerülnek forgalomba. A tálcára illeszkedő szőnyegek száma nem erre a csomagméretre, hanem egyetlen szőnyegre vonatkozik.

45. táblázat: Solaris 2000 tálcákhoz választható öntapadós szőnyegek

Solaris 4000 tálcákhoz

	Cikkszám	Unive	rzális	Kétlépcsős
		18 x 30	36 x 24	18 x 30
Nagy tapadású szőnyeg, 200 x 200 mm, csomag. egység: 2 db.	75004126	6	12	12
Nagy tapadású szőnyeg, 280 x 170 mm, csomag. egység: 2 db.	75004127	6	9	12
Kis tapadású szőnyeg, 200 x 200 mm, csomag. egység: 2 db.	75004111	6	12	12
Kis tapadású szőnyeg, 280 x 170 mm, csomag. egység: 2 db.	75004117	6	9	12

A szőnyegek 2 db-os kiszerelésben kerülnek forgalomba. A tálcára illeszkedő szőnyegek száma nem erre a csomagméretre, hanem egyetlen szőnyegre vonatkozik.

46. táblázat: Solaris 4000 tálcákhoz választható öntapadós szőnyegek

Tartozékok

Solaris 2000 I / 2000 R tálcákhoz

	Cikkszám	Univerzális
		11 x 14
Nagy tapadású szőnyeg, 200 x 200 mm, csomag. egység: 2 db.	75004126	1
Nagy tapadású szőnyeg, 280 x 170 mm, csomag. egység: 2 db.	75004127	2
Kis tapadású szőnyeg, 200 x 200 mm, csomag. egység: 2 db.	75004111	1
Kis tapadású szőnyeg, 280 x 170 mm, csomag. egység: 2 db.	75004117	2

A szőnyegek 2 db-os kiszerelésben kerülnek forgalomba. A tálcára illeszkedő szőnyegek száma nem erre a csomagméretre, hanem egyetlen szőnyegre vonatkozik.

47. táblázat: Solaris 2000 I/2000 R tálcákhoz választható öntapadós szőnyegek

Solaris 4000 I / 4000 R tálcákhoz

	Cikkszám	Univerzális
		18 x 18
Nagy tapadású szőnyeg, 200 x 200 mm, csomag. egység: 2 db.	75004126	4
Nagy tapadású szőnyeg, 280 x 170 mm, csomag. egység: 2 db.	75004127	2
Kis tapadású szőnyeg, 200 x 200 mm, csomag. egység: 2 db.	75004111	4
Kis tapadású szőnyeg, 280 x 170 mm, csomag. egység: 2 db.	75004117	2

A szőnyegek 2 db-os kiszerelésben kerülnek forgalomba. A tálcára illeszkedő szőnyegek száma nem erre a csomagméretre, hanem egyetlen szőnyegre vonatkozik.

48. táblázat: Solaris 4000 I/4000 R tálcákhoz választható öntapadós szőnyegek

Tartozékok

1. 2. 12. Általános tartozékok

1	1. 2. 12. Altalános tartozékok							
		Cikkszám						
	Felhordó tekercs öntapadós szőnyegekhez	75004124						
	Csavarhúzó a tartozékokhoz	75004131						
	Gázelosztó (Solaris 2000 I/2000 R)	SK2000-8GM						
	Gázelosztó (Solaris 4000 I/4000 R)	SK4000-8GM						
	Lombiktartó pótalkatrész készlet (csavarok)	SK0010						
	Pótalkatrészkészlet Solaris 2000, 2000 I/R és 4000 I/R tálcákhoz (tálcacsavarok, szerszám)	SK0100						
	Tálca pótalkatrész készlet SK4000-hez (tálca csavarok, szerszám)	SK0101						

49. táblázat: Választható általános tartozékok

1. 3. Szabványok és irányelvek

Régió	Irányelv	Szabványok
Európa	2006/42/EU	EN 61010-1
	Gépek irányelv	IEC 61010-2-051
	2011/65/EU RoHS	EN 61326-1, B osztály
	Bizonyos veszélyes anyagok	EN ISO 14971
	használatának korlátozása elektromos	EN ISO 9001
	Védelmi célok:	Kizárólag Solaris 2000 I/2000 R/
	2014/35/ELLKisfeszültségű irányely	4000 I/4000 R tálcákhoz:
		IEC 61010-2-010
	összeférhetőség (EMV)	
Észak-Amerika		ANSI/UL 61010-1
		IEC 61010-2-051
		EN 61326-1, B osztály
		CFR 47 FCC 15 EMV
		EN ISO 14971
		EN ISO 9001
		Kizárólag Solaris 2000 I/2000 R/
		4000 I/4000 R tálcákhoz:
		IEC 61010-2-010
Japán		IEC 61010-1
		IEC 61010-2-051
		EN 61326-1, B osztály
		EN ISO 14971
		EN ISO 9001
		<u>Kizárólag Solaris 2000 I/2000 R/</u> 4000 I/4000 R tálcákhoz: IEC 61010-2-010

50. táblázat: Szabványok és irányelvek

MEGJEGYZÉS: A készüléket bevizsgáltuk és megfelelt a digitális készülékek határértékeinek, az FCC előírások 15. része szerinti A osztályban. A határértékek megfelelő védelmet biztosítanak a káros zavarok ellen, ha a készüléket kereskedelmi környezetben használják. A készülék nagyfrekvenciás energiát állít elő és használ, amit ki is sugározhat, ha nem a kezelési útmutatójának megfelelően telepítik és használják, ezek a káros zavarok a rádiós kommunikációt érik. A berendezés lakóövezetben történő használata káros zavarokat eredményezhet. Ilyen esetben a felhasználó köteles saját költségére megszűntetni a zavarást.

1.4. Termék áttekintés

1. 4. 1. Solaris 2000



① Tálca szerelőlap; ② Grafikus felhasználói felület; ③ Hálózati kapcsoló; ④ Hálózati csatlakozó;
⑤ Biztosíték; ⑥ Ethernet-csatlakozó; ⑦ USB-csatlakozó

1. ábra: Solaris 2000 áttekintés

1. 4. 2. Solaris 4000



① Tálca szerelőlap; ② Grafikus felhasználói felület; ③ Hálózati kapcsoló; ④ Hálózati csatlakozó;
⑤ Biztosíték (visszaállítható); ⑥ Ethernet-csatlakozó; ⑦ USB-csatlakozó

2. ábra: Solaris 4000 áttekintés

Termék áttekintés

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



① Tálca szerelőlap; ② Grafikus felhasználói felület; ③ Hálózati kapcsoló; ④ Tömszelence; ⑤ Hálózati csatlakozó; ⑥ Biztosíték (visszaállítható); ⑦ Ethernet-csatlakozó; ⑧ USB-csatlakozó; ⑨ Burkolat

3. ábra: Solaris 2000 I / 2000 R áttekintés

1. 4. 4. Solaris 4000 I / 4000 R



① Tálca szerelőlap; ② Grafikus felhasználói felület; ③ Hálózati kapcsoló; ④ Tömszelence; ⑤ Hálózati csatlakozó; ⑥ Biztosíték (visszaállítható); ⑦ Ethernet-csatlakozó; ⑧ USB-csatlakozó; ⑨ Burkolat

4. ábra: Solaris 4000 I / 4000 R áttekintés

1. 4. 5. Csatlakozások

Nyitott rázógép



1) Hálózati csatlakozó; 2) Biztosíték; 3) Ethernet-csatlakozó

5. ábra: Hátulnézet (bal: Solaris 4000, jobbra: Solaris 2000)

Inkubált rázógép (csak fűtött) és hűtött rázógép (hűtött és fűtött)



1Hálózati csatlakozó; 2Biztosíték; 3Ethernet-csatlakozó; 4Leeresztő tömlő

6. ábra: Hátulnézet (bal: Solaris 2000 I / 2000 R, jobbra: Solaris 4000 I / 4000 R)

Hálózati csatlakozás

A rázógéphez olyan áramforrás szükséges, amely megfelel a specifikációinak. A hálózati kábel a csomagban található.

Cikkszám	Rázó	Műszaki adatok
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

7. ábra: A rázógépek tápellátására vonatkozó előírások

FIGYELMEZTETÉS A nem megfelelő hálózati csatlakozás vagy csatlakozó károsodást okoz. Ügyeljen arra, hogy a rázógépet csak olyan aljzatokhoz csatlakoztassa hozzá, amelyek a vonatkozó elektromos biztonsági szabványoknak megfelelő földeléssel rendelkeznek. Ne használja a rázót sérült vagy alulméretezett hálózati kábellel.

A hálózati csatlakozó mindig legyen hozzáférhető.

Az áramütés elkerülésére biztosítsa, hogy a rázó körüli terület száraz legyen. Folyadékok véletlen kiömlése vagy kifröccsenése esetén a folytatás előtt válassza le a rázógépet a hálózatról, távolítsa el a kiömlött folyadékot, és szüntessen meg mindenféle egészségügyi vagy biológiai veszélyt.

Ha nem használja a rázót, válassza le a hálózatról.

<u>Biztosíték</u>

A túláram következtében kioldott biztosíték visszaállítható. A rázó csak akkor kapcsolható be ismét, ha a biztosítékot visszaállították a megfelelő helyzetébe.

Ethernet

A rázón RJ-45 Ethernet csatlakozó van, amivel helyi hálózatra (LAN) kapcsolható. Kizárólag az IEC 60950-1 szabványnak megfelelő készülékek használhatók RJ45 Ethernet csatlakozóval. A később elérhető szoftverfrissítések az RJ45 Ethernet-csatlakozón keresztül telepíthetők.

<u>Tömszelencék</u>

A hőmérséklet-szabályozott rázógépek két oldalsó, tömszelencének nevezett nyílással rendelkeznek. Ezek a tömszelencék általában kupakkal vannak lezárva. A kupak levételét követően további eszközöket, például még egy hőmérsékletérzékelőt is üzembe helyezhet, vagy átvezetheti rajta a kábeleket vagy tömlőket, például a gázelosztó tömlőit.

<u>USB</u>

A rázógép két, A típusú USB 2.0 aljzatához szabványos USB-eszközök csatlakoztathatók. Kizárólag az IEC 60950-1 szabványnak megfelelő készülékek használhatók USB csatlakozóval.

Belső túlmelegedés elleni védelem

A hőmérséklet-szabályozott rázógépek a fűtőelemek mellé felszerelt, kézi visszaállítású termosztáttal is rendelkeznek. Ez a termikus eszköz kikapcsolja a fűtőelemeket, hogy megakadályozza a túlzott hőmérsékletet a rázógép burkolatán belül, ha meghibásodnának a keringető ventilátorok.

A kézi visszaállítású termosztátot kézzel kell visszaállítani a rázógép újbóli működése érdekében. Ezt kizárólag a Thermo Fisher Scientific ügyfélszolgálata hajthatja végre.

2. Szállítás és telepítés



Az üzemeltető felelős a rázó szabályos felállításáért.

MEGJEGYZÉS

Kiszállításkor azonnal ellenőrizze a csomagolás karton dobozát. Átvételkor azonnal ellenőrizze a szállítási sérüléseket, mielőtt a szállítmányt kicsomagolná. Ha sérülést észlel, rögzíttesse a szállítólevél másolatán és írassa alá a szállítóval.

Óvatosan nyissa ki a kartondobozt, és ellenőrizze az összes részegység ("51. táblázat: A szállítmány tartalma") hiánytalan meglétét, mielőtt kidobja a csomagolóanyagot. Ha a kicsomagolás után sérülést észlel, jelezze a speditőrnek és kérje a kár kivizsgálását.

Fontos! Ha nem az átvételt követő néhány napon belül kéri a kár kivizsgálását, a szállítmányozó mentesül a kártérítési kötelezettsége alól. Kérjen kár kivizsgálást.

2.1. Kicsomagolás

A kicsomagoláskor használja a darabjegyzéket annak megállapítására, hogy hiánytalanul érkezett-e. Ne dobja ki a csomagolóanyagot addig, amíg nem ellenőrizte az alkatrészek hiánytalan meglétét!

A szállítmány tartalma

Cikk	Mennyiség
Rázó	1
Hálózati csatlakozó kábel	1
Univerzális tálca	1*
Tálcacsavarok	
– Solaris 4000	4*
– egyéb kivitelek	3*
Kezelési útmutató (nyomtatott, angol)	1
Kezelési útmutató USB pendrive-on	1
Rögzítő szerszám (T-markolatú kulcs)	1
* A 2000 I, 2000 R, 4000 I és 4000 R rázógépeknél a tálca rázógépbe. Sem a tálca, sem a csavarok nem kaphatók k	a gyárilag be van építve a ülön alkatrészként.

51. táblázat: A szállítmány tartalma

Ha nincs minden rész a csomagban, forduljon a Thermo Fischer Scientifichez.

2.2. Telepítési hely

VIGYÁZAT Az UV sugárzásnak kitett műanyag alkatrészek csökkenő stabilitása károsan befolyásolhatja a védelmet. Ne tegye ki közvetlen napsütésnek vagy más UV forrásnak a rázót és a műanyagból készült tartozékait.

A rázót állítsa sík felületű asztalra vagy padra, ami elbírja a működő rázó és tartozékai ill. a minták súlyát. A rázót állítsa olyan dugaszoló aljzat közelébe, ami megfelel a típustábla követelményeinek. Hagyjon helyet a légáramlásnak, a tartozékoknak és a kényelmes kezelésnek a készülék körül.

A készülék felállítása során vegye figyelembe az alábbi követelményeket:

- Nyitott rázógép esetén a mozgóalkatrészekkel szembeni megfelelő távolság betartása, valamint a zúzódásos sérüléseknek és a szomszédos berendezések károsodásának az elkerülése érdekében hagyjon szabadon 8 cm-nyi (3 colnyi) helyet a tálca mind a négy oldalán. A nyitott rázógép tálcája a ház fölé nyúlhat.
- A rázógépek rezgéseket okoznak. Ne tároljon érzékeny készülékeket vagy veszélyes tárgyakat ill. anyagokat a biztonsági zónában.



8. ábra: Nyitott rázógépek beállítása 8 cm-es (3 colos) biztonsági távolsággal

A megfelelő szellőzés és az optimális teljesítmény garantálása érdekében a hőmérsékletszabályozott rázógépek esetében a következő távolsági követelményeket kell betartani:

- A Solaris 2000 I és 4000 I (csak fűtött) inkubált rázógépeknél a burkolat mind a négy oldalán 8 cm-es (3 colos) biztonsági távolságot kell tartani, lásd 9. ábra.
- A Solaris 2000 R és 4000 R hűtött rázógépek (hűtött és fűtött) a háznak mind a négy oldalán szabadon hagyott 8 cm-es (3 colos) távolsággal üzemeltethetők, ha a burkolat felett van szabad tér, amint az a 9. ábra bal oldalán látható.

 Ha a szabad tér a hűtött rázógép (hűtött és fűtött) Solaris 2000 R és 4000 R fölé kerül beépítésre, a burkolat körüli távolságot mind a négy oldalon 30 cm-re (12 colra) kell növelni, ahogyan az a 9. ábra jobb oldalán látható.



9. ábra: Az inkubált (csak fűtött) (balra) és a hűtött rázógépek (hűtött és fűtött) (jobbra), valamint a készülék fölött szabad térrel rendelkező hűtött rázógépek (hűtött és fűtött) beállítási távolsága a bal oldali ábrán látható.

FIGYELMEZTETÉS A rázás közben tartsa távol a veszélyes anyagokat ettől a zónától.

Az aljzat legyen:

- » legyen tiszta,
- » stabil, masszív, kemény és rezgésmentes legyen,
- » tegye lehetővé a rázó vízszintes felállítását,
- » viselje el a rázó súlyát.
- A rázót csak beltérben szabad használni.
- A rázót ne tegye ki hőhatásnak és erős napsütésnek.
- A felállítási hely legyen mindig jól szellőztetett.
- A hálózati csatlakozó mindig legyen hozzáférhető.

2.3. Szállítás

VIGYÁZAT A rázó leesése sérülést okozhat. A rázót mindig két oldalán fogva emelje meg. Sose emelje meg a rázót az előlapjánál vagy a rászerelt tálcánál fogva. Soha ne az átlátszó burkolat fogantyújánál fogva emelje fel a hőmérséklet-szabályozott rázógépet!

VIGYÁZAT A rázógép biztonságos működését a házon belül lecsapódó víz is befolyásolja. Ez akkor fordul elő, ha a rázógépet hűvösebb helyről melegebb, párásabb levegőjű helyre viszik, például egy raktárhelyiségből vagy a szállítókocsiból a laboratóriumi helyiségbe. A használatba vétel előtt hagyja legalább 2 órán át száradni a készüléket.

MEGJEGYZÉS Nyitott kivitelű rázógépek esetében a rázógép elmozdítása előtt mindig távolítsa el a tálcát, a rakományt és az összes tartozékot. Ha nem szereli le a tálcát, károsodhat a tálca szerelőlapja vagy a rázó mechanizmus. Hőmérséklet-szabályozott rázógépeknél az is elegendő, ha a rakományt eltávolítják a tálcáról.

MEGJEGYZÉS A rázó az ütésektől megsérülhet. A rázót álló helyzetben, lehetőség szerint az eredeti csomagolásában szállítsa.

A rázógép kezelése

A rázógép kezelésénél ügyeljen arra, hogy:

- a rázógépet mindkét oldalról, de soha ne elölről és hátulról emelje fel;
- annyi embert alkalmazzon a szállításához, amennyit a rázógép súlya megkövetel, de soha ne kevesebbet, mint két ember (további információért lásd a "Műszaki adatok" a 11. oldalon).





10. ábra: A rázó kézi szállítása

FIGYELMEZTETÉS A rázót mindig két oldalán fogva emelje meg. Soha ne emelje fel a rázógépet elölről vagy hátulról! A rázógépek nehezek (további információkért lásd a "Műszaki adatok" a 11. oldalon). A rázógép felemeléséhez és szállításához legalább 2 személy szükséges.

VIGYÁZAT Soha ne az átlátszó burkolat fogantyújánál fogja meg a hőmérséklet-szabályozott rázógépet annak áthelyezésekor! A fogantyú letörhet, a rázógép pedig a földre eshet, és súlyos személyi sérülés történhet.

2.4. Vízszintezés

A rázó nem rendelkezik vízszintbe állító szerkezettel. A rázó megfelelő felállításhoz vízszintezett aljzat szükséges.

MEGJEGYZÉS Ne helyezzen távtartó alátéteket vagy más lapos tárgyakat a rázógép lábai és a tartófelület közé a rázógép vízszintes helyzetbe állításához.

2.5. Hálózati csatlakozás

A rázógéphez olyan áramforrás szükséges, amely megfelel a specifikációinak. A hálózati kábel a csomagban található. További információért lásd a 7. ábra a 47. oldalon.

FIGYELMEZTETÉS A nem megfelelő hálózati csatlakozás vagy csatlakozó károsodást okoz. Biztosítsa, hogy a rázót csak megfelelően földelt dugaszoló aljzathoz csatlakoztassák. Ne használja a rázót sérült vagy alulméretezett hálózati kábellel.

MEGJEGYZÉS Az elektromágneses sugárzás interferenciát okozhat a kijelzőn. Ez nem károsítja a készüléket, nem korlátozza annak működését, és nem változtatja azt meg. Az elektromágneses sugárzás okozta interferencia elkerülése érdekében ne vigyen mobil eszközöket, például mobiltelefonokat a készülék közvetlen közelébe. Ne üzemeltesse a készüléket más, nagy elektromos teljesítményű készülékekkel közös áramkörön! Ne működtessen több készüléket közös hálózati elosztóról!

A rázó hálózatra kapcsolásához az alábbiak szerint járjon el:

- 1. Kapcsolja ki a hálózati kapcsolót a jobb oldalon.
- 2. Győződjön meg arról, hogy a kábel megfelel-e az országában érvényes biztonsági előírásoknak.
- Győződjön meg arról, hogy a centrifuga típustábláján levő feszültség és frekvencia adatok megegyeznek-e a hálózatéval.

A hálózati csatlakozó mindig legyen hozzáférhető.

Az áramütés elkerülésére biztosítsa, hogy a rázó körüli terület száraz legyen. Folyadékok véletlen kiömlése vagy kifröccsenése esetén válassza le a rázógépet a hálózatról, és a folytatás előtt távolítsa el a kiömlött folyadékot.

Ha nem használja a rázót, válassza le a hálózatról.

2.6. Alap beállítás

A készülék használata előtt végezze el az alábbi lépéseket:

 Kapcsolja be a készüléket. Megjelenik a Thermo Scientific cég emblémája. Nyomja meg az Start Setup gombot.



11. ábra: Alapbeállítás lekérése

2. Válassza ki a képernyő kívánt nyelvét. Nyomja meg a Tovább gombot.



12. ábra: Alap beállítások - Nyelv

 A készülék nevét szükség esetén a Készülék megnevezése párbeszédablakban adhatja meg. Nyomja meg a Tovább gombot.

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

13. ábra: Alap beállítások - Készülék megnevezés

4. A szövegmezőben szükség esetén megadhatja a várost és az országot. Ha beír három karaktert, a város és ország szövegmezőben megjelenik a megfelelő javaslatok listája. Szükség esetén kiválaszthat egy javaslatot, majd megérintheti a **Tovább** gombot.

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

14. ábra: Alap beállítások - Régió

5. Válassza ki a megfelelő dátum formátumot. Nyomja meg a Tovább gombot.

Set Date D	splay Format
MN	IM/DD/YYYY
O DD,	/MMM/YYYY
O YY	YY/MMM/DD
Back	Next

15. ábra: Alap beállítások - Dátum formátum

 A beállítókerék forgókapcsolóit elforgatva válassza ki az aktuális dátumot. Nyomja meg a Tovább gombot.



16. ábra: Alap beállítások - Dátum beállítás

Alap beállítás

 Válassza ki a kívánt időformátumot, és a beállítókerék forgókapcsolóit elforgatva állítsa be a pontos időt. Nyomja meg a **Tovább** gombot.



17. ábra: Alap beállítások - Óra beállítás

 Csak inkubált rázógépek (csak fűtött) és hűtött rázógépek (hűtött és fűtött) esetén: Válassza ki a kívánt hőmérsékleti mértékegységet. Nyomja meg a Tovább gombot.

Tempera	ture Units
	°C
0	۴F
Back	Next

18. ábra: Alapbeállítás – hőmérsékleti mértékegység kiválasztása

 Csak inkubált rázógépek (csak fűtött) és hűtött rázógépek (hűtött és fűtött) esetén: Kattintson a kék nyílra a túlmelegedési riasztás és a vészleállítás határértékeinek beállításához. Nyomja meg a Tovább gombot.



19. ábra: Alapbeállítás - a túl magas és túl alacsony hőmérséklet riasztási határértékeinek beállítása

- Ismételje meg az eljárást az alacsony hőmérséklet riasztási határértékei esetében is. A folytatáshoz érintse meg a **Tovább** gombot.
- Ha rendszergazda szintű hozzáférési kód beállítása nélkül szeretne továbblépni, hagyja bejelölve a Hozzáférés-ellenőrzés nélkül beállítást, és a folytatáshoz érintse meg a Tovább gombot.

Ha ekkor rendszergazda szintű hozzáférési kódot szeretne beállítani, érintse meg a Hozzáférés-ellenőrzéssel elemet.

	Access	s Mode
•	Open Mode Select to allow full access to	all functionality for all users.
0	Secure Mode Select to require users to ent changes to stored parameter (All users still able to use prog	ter the admin passcode to make rs, programs, and settings. rams and set new run parameters.
1	Back	Next

20. ábra: Alapbeállítás - hozzáférési mód kiválasztása

 A hozzáférési kód kérése üzenet megjelenésekor írja be a rendszergazda szintű hozzáférési kódot a billentyűzet segítségével, majd érintse meg a **Tovább** gombot.



- 21. ábra: Alapbeállítás hozzáférési mód kiválasztása
- A hozzáférési kód kérése üzenet következő megjelenésekor adja meg a rendszergazda szintű hozzáférési kódot.
- A hozzáférési kód kérése üzenet harmadik megjelenésekor a megerősítéshez adja meg ismét a rendszergazda hozzáférési kódot.

		Adm	in Pass	code				Adm	in Pass	code	
f you do r	iot wis fields	h to change below blanl	the pa and ta	sscod p the	e at this Next bu	time, leave the tton.		Please re-ent	er the n	ew passcode.	
	*	*	*		*	*	*	*	*	*	*
		7	8	9				7	8	9	
		4	5	6				4	5	6	
		1	2	3				1	2	3	
		0	Cle	ear				0	Cle	ar	

22. ábra: Alapbeállítás - új rendszergazda szintű hozzáférési kód megadása és megerősítése

57 / 140

15. A megjelenő megerősítő képernyőn a folytatáshoz érintse meg a Tovább gombot.



23. ábra: Alapbeállítás – a rendszergazda szintű hozzáférési kód módosításának megerősítése

 Megjelenik egy képernyő a rázógép általános fizikai telepítési utasításaival. Nyomja meg a Tovább gombot.

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			

24. ábra: Alap beállítás - Telepítés és csatlakoztatás

 Megjelenik A konfigurálás befejeződött képernyő. A folyamat lezárásához érintse meg a Tovább gombot.

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

25. ábra: Alapbeállítás – Kalibrálás vagy az alapbeállítás befejezése

MEGJEGYZÉS Inkubált rázógépek (csak fűtött) és hűtött rázógépek (hűtött és fűtött) esetén a Kalibrálás gomb megérintésével hajthatja végre a folyamatot, amint azt a "Hőmérsékletkalibrálás" a 126. oldalon.

2.7. Tárolás



VIGYÁZAT

A rázó és tartozékainak tárolása előtt a teljes rendszert meg kell tisztítani, és ha szükséges fertőtleníteni ill. dekontaminálni. Ne helyezze raktárba a rázót és tartozékait mindaddig, amíg nem tisztázták egyértelműen a kontaminációs állapotát. Ha nem biztos az eljárással kapcsolatban, forduljon a Thermo Fisher Scientific ügyfélszolgálatához ("Tisztítás" a 123. oldalon, "Fertőtlenítés" a 124. oldalon és "Dekontaminálás" a 124. oldalon).

- A raktárba helyezés előtt meg kell tisztítani, és szükség esetén fertőtleníteni ill. dekontaminálni a rázót és tartozékait.
- Teljesen szárítsa meg a rázót és tartozékait, mielőtt raktárba helyezné.
- A rázót tiszta, pormentes helyen tárolja.
- A rázót a saját lábain állva tárolja.
- Ne tárolja a rázót közvetlen napsütésben.

2.8. Elküldés



VIGYÁZAT

A rázó és tartozékainak elküldése előtt a teljes rendszert meg kell tisztítani, és ha szükséges fertőtleníteni ill. dekontaminálni. Ne helyezze raktárba a rázót és tartozékait mindaddig, amíg nem tisztázták egyértelműen a kontaminációs állapotát. Ha nem biztos az eljárással kapcsolatban, forduljon a Thermo Fisher Scientific ügyfélszolgálatához ("Tisztítás" a 123. oldalon, "Fertőtlenítés" a 124. oldalon és "Dekontaminálás" a 124. oldalon).

A rázó elküldése előtt ügyeljen az alábbiakra:

- A rázót meg kell tisztítani ill. dekontaminálni.
- A dekontaminálást tanúsítvánnyal kell igazolni. A Thermo Fisher Scientific ügyfélszolgálatánál kérhet dekontaminációs tanúsítványt.

3. Kezelés

3.1. Be-/Kikapcsolás

A jobb oldalon levő hálózati kapcsolót nyomja meg a rázó be- vagy kikapcsolásához (I vagy 0).

Rendszerindulás közben a kezelőfelületen a Thermo Scientific logója látható.

Ha készen áll, a felhasználói felület megjeleníti a rázógép aktuális üzemállapotát.

3. 2. Grafikus felhasználói felület

A kezdőképernyő a rázógép grafikus felhasználói felületének alapértelmezett képernyője. Ez a rázógép üzemeltetésének kiindulási pontja.

Lehetőségek a kezdőképernyőn:

- a legfontosabb üzemi paramétereknek, például a rázás fordulatszámának, időtartamának és hőmérsékletének a beállítása; (kizárólag hőmérséklet-szabályozott rázógépek esetén)
- a rázógép elindítása és leállítása;
- az üzemállapotra vonatkozó információk megjelenítése, valamint a riasztási és figyelmeztető üzenetek szerkesztése;
- átváltás az üzemállapotra vonatkozó információkat és beállítási lehetőségeket megjelenítő egyéb képernyőkre.

A nyitott és a hőmérséklet-szabályozott rázógépek képernyőtartalma némileg eltérő.

A 26. ábran egy nyitott rázógép kezdőképernyője látható.



 Navigációs sor 2 Névleges fordulatszám 3 Start vagy Stop (4) Információ Időzítő - folyamatos üzem (5) vagy előre beállított időtartamú üzemmód (6) Tényleges fordulatszám Üzemállapot és statisztikai $\overline{(7)}$ adatok (8) Pontos idő (9) Készülék megnevezés

26. ábra: GUI Nyitott rázógép kezdőképernyője

A hőmérséklet-szabályozott rázógépek kezdőképernyőjén további hőmérsékletszabályozók találhatók, amint az a 27. ábran látható.



27. ábra: GUI Hőmérséklet-szabályozott rázógép kezdőképernyője

3. 2. 1. A rázógép legfontosabb üzemi paramétereinek beállítása

A kezdőképernyőn egyetlen érintéssel megjeleníthető a rázógép legfontosabb üzemi paramétereinek beállítási képernyője.

A fordulatszám beállítása



28. ábra: Névleges fordulatszám

 A kívánt fordulatszám beállításához érintse meg az egyes beállítókerekek tárcsái felett vagy alatt látható nyilakat. Vagy módosíthatja is a kívánt névleges fordulatszámot a beállítókerék egyes forgókapcsolóinak elforgatásával. 3. Nyomja meg a **Mentés** gombot.

MEGJEGYZÉS Ha olyan felugró ablak jelenik meg, amely arról tájékoztatja, hogy a névleges fordulatszám a megengedett tartományon kívül van, akkor az Ön beállítása a rázógép által támogatott üzemi fordulatszám-tartományon kívül van, amint azt a "Műszaki adatok" a 11. oldalon olvashatja. Helyesbítse a beállítást, és folytassa.

 Nyugtázza az ezután megjelenő Megtörtént a beállításainak mentése párbeszédablakot a főképernyőre történő visszatéréshez.

Körpálya számítógép: a hozzávetőleges fordulatszám kiszámítása a készülék előző beállításai alapján

A Névleges fordulatszám párbeszédablak mezőben a Fordulatszám kiszámítása elem megérintésével is megnyithatja a Körpálya számítógép képernyőt: A körpálya számítógép akkor hasznos, ha egy eltérő mechanikai felépítésű rázógépről átvált a jelen kezelési útmutatóban leírt Solaris sorozatú rázógépek valamelyikére. Ez a számítógép a körpálya méretének különbségét használja fel ahhoz, hogy nagyjából megbecsülje a hasonló eredmény eléréséhez szükséges névleges fordulatszámot.

MEGJEGYZÉS Ne használja fel azonnal a körpálya számítógéppel kapott eredményt a gyártási mintákhoz, hanem végezzen el egy sor tesztet megerősítés céljából! A sejtek gyorsabban vagy lassabban növekedhetnek, más fehérjéket képezhetnek, vagy megsérülhetnek (vagy elpusztulhatnak) a jelenlévő nyíróerőktől függően, mivel a fizikai feltételek megváltoznak egy másik körpályára való mozgás közben.



1. Válasszon mértékegységet (mm vagy Zoll).

29. ábra: Körpálya számítógép

 Válassza ki az előző körpálya nagyságát. Nyomja meg a Számolás gombot. A fordulatszám kiszámításához az alábbi ablak jelenik meg.



30. ábra: Körpálya kiszámítva

- 3. Nyomja meg az **OK** gombot.
- Kiszámíthatja az ajánlott fordulatszámot egy vevőspecifikus körpálya mérethez. Ügyfélspecifikus körpályára vonatkozó számítás elvégzéséhez válassza az Egyéb lehetőséget.
- 5. Adja meg a körpálya méretét.



31. ábra: Ügyfélspecifikus körpálya méretének kiszámítása

6. Nyomja meg a Számolás gombot.

Megjelenik a Fordulatszám kiszámítása ablak és a megadott nagysághoz tartozó fordulatszámot mutatja.

Működési idő beállítása

A rázógép folyamatos üzemmódban vagy előre beállított időtartammal üzemeltethető. Folyamatos üzemben a rázógépet szükség esetén kézzel is leállíthatja. Előre beállított időtartamú üzemmódban a rázógép hajtása automatikusan leáll, amikor lejár az időzítő. Az időzítőt a következőképpen jelenítheti meg:

- eltelt idő: ennyi ideje fut a rázógép a Start gomb megnyomása óta, vagy
- hátralévő idő: ennyi ideig fut még a rázógép, amíg le nem jár az időzítő

- Érintse meg az Idő mezőt (⑤ a 26. ábran és 27. ábran) a 32. ábran látható Időmód képernyő megnyitásához.
- 2. Válassza az Üzemelés előre beállított időtartammal módot.



32. ábra: Időmód kiválasztása

- 3. Az Idő beállításnál adja meg az órát és percet.
- 4. Nyomja meg a Mentés gombot.
- A főképernyőre történő visszatéréshez érintse meg a megjelenő kérést, ami a Megtörtént a beállításainak mentése nyugtázására szolgál.

A hőmérséklet beállítása

Hőmérséklet-szabályozott rázógépek esetén a Hőmérséklet névleges értéke mezővel bármikor beállíthatja az alkalmazáshoz szükséges hőmérsékletet.

- 2. Érintse meg a beállítókerék egyes forgókapcsolói feletti vagy alatti nyilakat annak az alapértelmezett hőmérsékletnek az előzetes beállításához, amelyet a rázógépnek fenn kell tartania. A kívánt hőmérsékletet a beállítókerék egyes forgókapcsolóinak külön-külön történő elforgatásával is megváltoztathatja.



33. ábra: A hőmérséklet névleges értékének beállítása: Képernyő kalibrációs kijelzővel

3. Nyomja meg a **Mentés** gombot.

MEGJEGYZÉS Ha egy felugró ablak jelenik meg, amely arról tájékoztatja, hogy a hőmérséklet névleges értéke a megengedett tartományon kívül van, akkor az Ön beállítása kívül esik a rázógép által támogatott üzemi hőmérséklettartományon, amint azt a "Műszaki adatok" a 11. oldalon jelzi. Helyesbítse a beállítást, és folytassa.

MEGJEGYZÉS Ha egy felugró ablak jelenik meg, amely arról tájékoztatja, hogy a kiválasztott hőmérséklet az aktuális környezeti hőmérséklet következtében a megengedett tartományon kívül esik, akkor a környezeti hőmérséklet kívül esik a rázógép üzemi hőmérséklettartományán, ahogyan az a "Műszaki adatok" a 11. oldalon szerepel. Érintse meg az **OK** gombot az értesítés nyugtázásához és a folytatáshoz, vagy válasszon másik hőmérsékletet.

MEGJEGYZÉS A Hőmérséklet névleges értéke képernyőn megjelenik a kalibrálát hőmérséklet és a kalibrálás során beállított eltolás, amint az a 33. ábra bal alsó részén látható. A kalibrálási folyamatot a "Hőmérsékletkalibrálás" a 126. oldalon ismertetjük.

 A főképernyőre történő visszatéréshez érintse meg a megjelenő kérést, ami a Megtörtént a beállításainak mentése nyugtázására szolgál.

MEGJEGYZÉS A hőmérséklet-szabályozott rázógépek a beállított hőmérséklet elérésekor meghatározott módon viselkednek, amit "túllengésnek" nevezünk. A hőmérséklet túllengés azt jelenti, hogy a kamra névleges hőmérsékletét először kissé magasabb (vagy alacsonyabb) hőmérsékletre vezérli a rendszer, majd megközelíti a névleges hőmérsékletet. A névleges hőmérséklethez közeledve a kijelző nem jeleníti meg ezt a túllengést. Ehelyett a kamrahőmérséklet növekedése (vagy csökkenése) jelenik meg a névleges hőmérséklet eléréséig.

A rázógép elindítása és leállítása

1. A rázógép elindításához érintse meg a Start gombot.

A Start gomb Szünet gombra vált.



34. ábra: Start és Stop gomb

2. A rázógép leállításához érintse meg a Stop gombot.

3. 2. 2. Üzemállapot

Ha a rázógép üzemállapota megfelelő, a kezelőmezőn, az üzemállapot és a statisztikai adatok



területen (⑦ a 26. ábran és ⑧ a 27. ábran) egy zöld szív szimbólum jelenik meg. A zöld szív szimbólum megérintésekor megnyílik az Üzemállapot képernyő. Az Üzemállapot képernyő statisztikai adatokat tartalmaz a rázógép működéséről, a rázási és üzemórák

összesített számát, illetve az utolsó munkamenetet is beleértve. A hőmérséklet-szabályozott rázógépek a hűtési és fűtési folyamatok óraszámait is kijelzik.



35. ábra: A rázógép statisztikai adatai: Nyitott rázógép (balra) és hőmérséklet-szabályozott rázógép (jobbra)

Figyelmeztetés

Ha figyelmeztető üzenet jelenik meg, a kezelőmező egy gördülő szöveggel rendelkező, sárga sávot

jelenít meg az aktuális képernyő felső részén. Ezenkívül egy riasztási hangjelzés is megszólal. A sárga figyelmeztető sáv eltűnik a gördülő szöveggel együtt, miután az üzenet kétszer lefutott. Egyedül a sárga figyelmeztető háromszög jelzi azt, hogy egy



vagy több riasztási üzenet van-e a rázógépre vonatkozóan. A figyelmeztető háromszögen belüli kék kör fehér szegéllyel azt jelzi, hogy hány aktív figyelmeztetés áll fenn. Az Információ és üzemállapot területen található figyelmeztető háromszög (⑦ a 26. ábran és ⑧ a 27. ábran) megérintésével megnyithatja az összes, éppen aktív figyelmeztető üzeneteket tartalmazó képernyőt. A legutóbbi figyelmeztetés részletesen jelenik meg, így az összes információt láthatja, amint az a 36. ábran látható. A listát görgetheti és bármelyik bejegyzést megérintve megnyithatja annak részletes formáját.



36. ábra: Figyelmeztető üzenetek listája

Az aktív figyelmeztető üzenetet a figyelmeztető üzenet bejegyzés melletti jelölőnégyzet megérintésével választhatja ki. A **Nyugtázás** gomb megérintése után a rendszer megkísérli törölni a figyelmeztető üzenetet a listából. Ha az összes figyelmeztető üzenet törlődött, az üzemállapot szimbólum visszaváltozik az eredeti zöld szívre.

<u>Riasztás</u>

Ha a készülék riasztást ad ki, a rázó azonnal leáll, hogy elkerülje a minták és/vagy magának a



készüléknek a sérülését. A riasztási üzenetet nyugtáznia kell a kezelőmezőben ahhoz, hogy folytathassa az üzemeltetést.

Ha riasztást ad ki, a kezelőfelület pillanatnyi képernyőjének tetején piros sáv jelenik meg. Az Információ és üzemállapot területen (⑦ a 26. ábran és ⑧ a 27. ábran) egy piros riasztási csengő jelenik meg, mindkét oldalán hanghullámokkal. Megszólal egy folyamatos riasztó hang is.

A piros riasztási sáv alatt egy görgethető szöveges üzenet jeleníti meg az aktuális riasztási üzenet összefoglalóját. A jobb oldalon megjelenik egy szundi gomb, amellyel átmenetileg lenémíthatja a riasztási üzenetet. Ha a némítási idő alatt nem hárítják el a riasztás okát, ismét megszólal a hangjelzés. A némítás ideje a Beállításoknál adható meg. További információkért lásd a "Riasztások/ Figyelmeztetések" a 69. oldalon.

Miután megérintette a csengő ikont az Információ és üzemállapot területen (⑦ a 26. ábran és ⑧ a 27. ábran), megjelenik egy képernyő, amely az éppen aktív riasztási üzenet összes részletét megmutatja, amint az a 37. ábran látható. A listát görgetheti és bármelyik bejegyzést megérintve megnyithatja annak részletes formáját.



37. ábra: Riasztási üzenetek listája

Az aktív riasztási üzenetet a riasztási bejegyzés melletti jelölőnégyzet megérintésével választhatja ki. A **Nyugtázás** gomb megérintésekor a rendszer megkísérli törölni a riasztási üzenetet a listából, ha megszüntette az okot. Amikor törlődött az összes riasztási üzenet, az üzemállapot szimbólum visszaváltozik az eredeti zöld szívre.

<u>Hiba</u>

Meghibásodás esetén a rázógép hibaüzenetet jelenít meg, és azonnal leáll, hogy elkerülje a minták és/vagy magának a készüléknek a károsodását. A képernyő pirosra vált, és nincs lehetőség további beavatkozásra. Hibaüzenet jelenik meg egy hibakóddal együtt, ahogy a 38. ábra példája mutatja.



38. ábra: Hibajelző képernyő

Javítsa ki a problémát az alábbiak szerint, és kísérelje meg az üzemeltetés folytatását:

- 1. Jegyezze fel a hibajelző képernyőn látható hibakódot.
- 2. A riasztási hang némításához érintse meg a Nyugtázás gombot.
- 3. Ki- és visszakapcsolással indítsa újra a rázógépet.
- Ha a hibaüzenet továbbra is látható, lépjen kapcsolatba a Thermo Fisher Scientific ügyfélszolgálattal, és tájékoztassa őket a hibajelző képernyőn megjelenő hibakódról.

3. 2. 3. Beállítások

A navigációs sáv második ikonja a beállításoké. A **Beállítások** elem megérintésekor az alábbi képernyő jelenik meg. A Beállítások képernyő több gombot tartalmaz, mint amennyi egyszerre megjeleníthető a kijelzőn. A többi gomb megjelenítéséhez le kell görgetnie, ahogyan az a 39. ábra jobb oldalán látható.





39. ábra: Képernyőbeállítások

Riasztások/Figyelmeztetések

A Riasztások/Figyelmeztetések képernyőn meghatározhatja, hogy a riasztási és figyelmeztető üzenetek hogyan és mikor jelenjenek meg.



40. ábra: A riasztási és figyelmeztető üzenetek előzetes beállításainak megadása

Hangerő: Ezzel a csúszkával a riasztás vagy figyelmeztetés esetén kiadott hang erősségét módosíthatja. Húzza a csúszkát balra a hangerő csökkentéséhez, vagy jobbra a hangerő növeléséhez, majd érintse meg a **Mentés** elemet. Ez a hangerő-beállítás a riasztási és a figyelmeztető üzeneteket is érinti.



41. ábra: A riasztási és figyelmeztető üzenetek hangerejének beállítása

Túl magas hőmérséklet/túl alacsony hőmérséklet riasztás (kizárólag hőmérsékletszabályozott rázógépeknél): Ezzel a két beállítással a hőmérséklet meghatározott névleges értékéhez viszonyítva állíthatja be a túl magas és túl alacsony hőmérséklet riasztási határértékét (lásd "A hőmérséklet beállítása" a 64. oldalon). Amint a rázókamrában a hőmérséklet a túl alacsony hőmérséklet határértéke alá csökken, vagy túllépi a túl magas hőmérséklet határértékét, a rázógép riasztási üzenetet jelenít meg. Válasszon hőmérséklet határértéket, és érintse meg a Mentés elemet.



42. ábra: Riasztási határértékek beállítása túl magas és túl alacsony hőmérséklethez

MEGJEGYZÉS A riasztási határértékek ennek megfelelően változnak, ha módosítja a hőmérséklet névleges értékét. Lekapcsolás túl magas hőmérséklet/túl alacsony hőmérséklet esetén (kizárólag hőmérséklet szabályozott rázógépek esetén): Ezzel a két beállítással a hőmérséklet meghatározott névleges értékéhez viszonyítva állíthatja be a biztonsági lekapcsolással kapcsolatos túl magas és túl alacsony hőmérséklet határértékét (lásd "A hőmérséklet beállítása" a 64. oldalon). Amint a rázókamrában a hőmérséklet a túl alacsony hőmérséklet határértéke alá csökken, vagy a túl magas hőmérséklet határértéke fölé emelkedik, a rázógép automatikusan lekapcsol a kamrában feldolgozott minták védelme érdekében. Válasszon ki egy határértéket, és érintse meg a **Mentés** gombot.

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

43. ábra: Biztonsági lekapcsolással kapcsolatos túl magas és túl alacsony hőmérséklet határértékek beállítása

MEGJEGYZÉS A hiba határértékek ennek megfelelően eltolódnak, ha módosítja a hőmérséklet névleges értékét.

Némítás időtartama: A némítás időtartama határozza meg azt, hogy mennyi időre némuljon el egy riasztási üzenet akkor, ha megnyomják a főablak felső részén található piros sávban lévő Szundi gombot (lásd "Riasztás" a 67. oldalon). A némítási időt 5, 10 vagy 15 perce állíthatja be. Az alapbeállítás 10 perc. A módosítások mentéséhez érintse meg a **Mentés** gombot.



44. ábra: A riasztási üzenetek némítási időtartamának beállítása

Kalibrációs üzenetek letiltása (kizárólag hőmérséklet-szabályozott rázógépek esetén): Ezekkel a beállításokkal letilthatja a rázógépén a hőmérsékletmérés kalibrálására rendszeresen emlékeztető üzeneteket (további információkért lásd a "Hőmérsékletkalibrálás" a 126. oldalon). Kiválaszthatja, hogy a kalibrálási emlékeztetőket teljesen vagy csak egy program végrehajtása közben kívánja-e letiltani.



Hozzáférés ellenőrzés

A Hozzáférés-ellenőrzés képernyőn rendszergazda szintű hozzáférési kódot rendelhet hozzá a korlátozott hozzáférési jogokkal rendelkező folyamatokba történő belépéshez.

A rázógép alapkivitel szerint úgy van beállítva, hogy hozzáférés-ellenőrzés nélkül működjön, azaz a készülékhez történő hozzáféréshez és működtetéshez nem szükséges belépési kódot megadni. A "Hozzáférés-ellenőrzéssel" üzemmódra történő átváltáshoz minden olyan felhasználónak szüksége van a rendszergazda szintű hozzáférési kódra, aki módosítani kívánja a rázógép beállításait. A Hozzáférés ellenőrzése opció választása esetén a kijelző jobb felső sarkában Bejelentkezés gomb jelenik meg.

Az alábbi összehasonlítás mutatja, hogy mikor van szükség a hozzáférési kódra a működtetéshez.

Teendő	Hozzáférési kódra van szükség a hozzáférés ellenőrzése nélküli módban	A hozzáférést ellenőrző módban hozzáférési kód szükséges
A fordulatszám, az időtartam és a hőmérséklet névleges értékeinek beállítása	Nem	Nem
A rázógép elindítása és lefékezése	Nem	Nem
Körpályák kiszámítása	Nem	Nem
Riasztási és figyelmeztető üzenetek nyugtázása	Nem	Nem
Képernyőfájlok és információk megjelenítése	Nem	Nem
Üzemállapot kijelzése	Nem	Nem
Eseménynapló és diagramok megtekintése és exportálása	Nem	Nem
Riasztások átmeneti némítása	Nem	Nem
Programok használata	Nem	Nem
Programok létrehozása, szerkesztése vagy törlése	Nem	lgen
Kijelző beállítások módosítása	Nem	lgen
Konfigurációs beállítások módosítása	Nem	lgen
Riasztási és figyelmeztető beállítások módosítása	Nem	lgen
A hozzáférés-ellenőrzés beállításainak módosítása	Nem	lgen
Programok importálása vagy exportálása	Nem	lgen
Gyári beállítások visszaállítása	Nem	lgen
Kalibrálás végrehajtása	Nem	lgen
Készülékszoftver-frissítés telepítése	lgen	lgen

52. táblázat: Hozzáférési kód kötelező megadása ellenőrzött vagy ellenőrizetlen módban

A 46. ábran a Hozzáférés-ellenőrzés képernyő látható hozzáférés-ellenőrzés nélkül és hozzáférésellenőrzéssel.



46. ábra: Hozzáférés-ellenőrzés: Hozzáférés-ellenőrzés nélkül (balra) és hozzáférés-ellenőrzéssel (jobbra)

Ha megérinti a **Hozzáférés-ellenőrzéssel**, majd pedig a **Mentés** elemet, a módosítás megerősítése előtt a rendszer kérni fogja a rendszergazda szintű hozzáférési kódot. A gyárban az összes készülékbe azonos adminisztrátori kódot programozunk. Ez a kód a kézikönyvben található.

Hasonlóképpen, bármely felhasználó, aki hozzáférési kóddal védett beállítást kísérel megváltoztatni, a rendszergazda szintű hozzáférési kódot kell hogy megadja.

MEGJEGYZÉS Az adminisztrátori kód gyárilag 00000 beállítású.

A Rendszergazda szintű hozzáférési kód módosítása gombbal módosíthatja az előre beállított hozzáférési kódot. A Rendszergazda szintű hozzáférési kód módosítása gomb megérintésekor a rendszer megkéri, hogy előbb adja meg az aktuális hozzáférési kódot, majd az új hozzáférési kódot, azután pedig megerősítés céljából ismét az új hozzáférési kódot. Az Új hozzáférési kódjának mentése megtörtént üzenet azt jelzi, hogy a művelet sikeres volt.

<u>Kalibrálás</u>

A Kalibrálás képernyő (kizárólag hőmérséklet-szabályozott rázógépek esetén) egy sor olyan képernyőt nyit meg, amelyek lehetővé teszik a rázógép kalibrálását a pontos hőmérsékletszabályozás érdekében. Ezt a műveletet a "Hőmérsékletkalibrálás" a 126. oldalon.
<u>Üzemi paraméterek</u>

Az Üzemi paraméterek képernyő lehetővé teszi azoknak az alapértelmezett üzemi paramétereknek az előzetes beállítását, amelyekkel a készüléket el kell indítani:

Névleges fordulatszám: A fordulatszám névleges értékét 15 és 525 fordulat/perc közötti értékre állíthatja be. A fordulatszám beállításához forgassa el a beállítókerék forgókapcsolóját, majd érintse meg a Mentés gombot. A fordulatszám beállítási képernyő használatára vonatkozó részletes utasításokat lásd a "A fordulatszám beállítása" a 61. oldalon.

Időmód: A rázógép folyamatos üzemmódban vagy előre beállított időtartammal üzemeltethető. Folyamatos üzemben a rázógépet szükség esetén kézzel is leállíthatja. Előre beállított időtartamú üzemmódban a rázógép hajtása automatikusan leáll, amikor lejár az időzítő. Előre beállított időtartamú működés esetén előre beállíthat egy alapértelmezett futási időt, és kiválaszthatja, hogy a felhasználói felület az eltelt időt vagy a hátralévő időt jelenítse-e meg a rázógép működése közben. Az Időmód képernyő használatára vonatkozó részletes utasításokat lásd az "Működési idő beállítása" a 63. oldalon.

Hőmérséklet névleges értéke (kizárólag hőmérséklet-szabályozott rázógépek esetén): Ez a beállítás lehetővé teszi, hogy beállítson egy olyan alapértelmezett hőmérsékletet, amelyet a rázógépnek indításkor meg kell jelenítenie. A Hőmérséklet névleges értéke képernyő használatára vonatkozó részletes utasításokat lásd a "A hőmérséklet beállítása" a 64. oldalon.

Automatikus újraindítás: Ezzel a funkcióval újraindíthatja a készüléket normál elindítás/leállítás, program végrehajtása vagy hőmérsékletkalibrálás közben bekövetkezett áramkimaradás után. Ha az Automatikus újraindítás opció beállítása **Nem**, áramkimaradás után a készülék nem indul újra.

<u>Kijelző</u>

A Kijelző beállításoknál tudja a különböző megjelenítési lehetőségeket beállítani.

Fényerő: A kijelző fényerejének beállításához használja a csúszkát vagy a + / - gombokat.



47. ábra: A képernyő fényerejének beállítása

Nyelv: A kijelző nyelvének megváltoztatásához érintse meg a **Nyelv** gombot. Válassza ki a kívánt nyelvet a beállítókerék elforgatásával, majd érintse meg a **Mentés** gombot.



48. ábra: A kijelző nyelvének kiválasztása

Mértékegységek (kizárólag hőmérséklet-szabályozott rázógépek esetén): érintse meg a Mértékegységek gombot, hogy átváltsa a hőmérséklet-kijelzéseket Celsius fok °C és Fahrenheit °F között a grafikus felhasználói felület képernyőin.

Units of Measure	°C 🔨
	°F

49. ábra: A hőmérséklet-kijelzés mértékegységének kiválasztása

Dátum: A dátum beállításához érintse meg a Dátum gombot. Forgassa el a beállítókerék forgókapcsolóit a hónap, a nap és az év kiválasztásához, majd érintse meg a **Mentés** gombot.



50. ábra: Hónap, dátum és idő beállítása

Dátum formátum: A dátumformátum beállításához érintse meg a Dátumformátum gombot. Érintse meg a kiválasztásnak megfelelő dátumformátum kiválasztó jelölőnégyzetét (például: HH/NN/ÉÉÉÉ megjelenítése, mint 2015. április 15.), majd érintse meg a Mentés gombot.



51. ábra: Dátumformátum beállítása

Idő: Az idő és az időformátum beállításához érintse meg az Idő gombot. Érintse meg a 12 óra vagy a 24 óra gombot, forgassa el a beállítókerék óra, perc és AM/FM (csak 12 órás formátumban) forgókapcsolóit, majd érintse meg a Mentés gombot.



52. ábra: Az idő beállítása és kijelzése 12 órás vagy 24 órás óra formátumban

Alvó mód: Érintse meg az Alvó üzemmód elemet, hogy a rázógép kijelzője 15 perc inaktivitás után alvó üzemmódba lépjen. Alvó üzemmódban a képernyő az Aktiváláshoz érintse meg felhasználónak szóló felszólítást jeleníti meg.



53. ábra: Alvó üzemmód aktiválása

Régió: A készülék működési területének beállításához érintse meg a **Régió** gombot. Érintse meg a Város, ország szövegmezőt, és kezdje el beírni a város nevének első betűit. A felhasználói felület az első három karakter beírása után lehetséges találatokat javasol. Fogadja el a megjelenő lehetőségek egyikét, vagy folytassa a teljes név beírását, majd érintse meg a képernyőn megjelenő billentyűzeten a **Mentés** elemet.

< ♠	Sh 11 Re	ake La 22:24/ gion	b 1 AM	Enter the		2	Joseet to	where the	8 7000		C Shake Lab 1 11:22:24AM Region		
⇔ ₀ <u>₿</u>	Cit	y, Coui	ntry:	N		Journey c	ioscut to	your un			City, Country	Enter the city and country closest to your time zone New York, United States	
Q	w	E	R	T	Y	U	1	0	Ρ		Ē		
A	S	D	F	G	Н	J	К	L	S	ave			
1	z	x	С	V	В	N	м			1	-		
?!1	23								+			Save	()

54. ábra: Régió beállítása

Készülék megnevezés: A rázógép elnevezéséhez vagy átnevezéséhez érintse meg a Készülék megnevezése gombot. Érintse meg a Készülék megnevezése szövegmezőt, és kezdje el a gépelést. Ha végzett, érintse meg a kijelzőn megjelenő billentyűzeten a Mentés elemet.

< ♠	Sh 11 Ur	ake La :22:24) hit Nam	b 1 AM ie		(2					< ♠	Shake Lab 1 11:22:24AM Unit Name	\bigcirc	
٥.	U	nit Na	me	Sh	ake La	ab 1]	۰.			
Q	W	E	R	Т	Y	U	T	0	Ρ	•	Ë	Unit Name:	Shake Lab 1	
А	S	D	F	G	Н	J	К	L	Sa	ive				
1	Z	Х	С	۷	В	Ν	М	@		1	÷		_	
?!1	23							-	-	.			Save	()

55. ábra: Készülékmegnevezés hozzárendelése a rázógéphez

Ügyfélspecifikus testre szabás menü: A fő navigációs sáv alsó két ikonjának testre szabásához érintse meg az Ügyfélspecifikus testre szabás menü gombot. Húzza a kívánt szimbólumot a főképernyő területéről a lecserélni kívánt szimbólumra. A megerősítéshez érintse meg a Mentés gombot. MEGJEGYZÉS A Menü visszaállítása gombot megérintve bármikor visszaállíthatja a navigációs sávot a gyári beállításokra.



56. ábra: A navigációs sáv testre szabása

Fájlok és info

Ez a képernyő jeleníti meg a sorozatszámot, a GUI-alkalmazás szoftverét és annak verzióját, a rázógép fő vezérlő készülékszoftverét és annak verzióját, a paraméterfájlt és annak verzióját, valamint az aktuális operációs rendszert.

A **Gyári beállítások visszaállítása** elem megérintésével visszaállíthatja a rázógép gyári beállításait. A gyári beállítások visszaállításához rendszergazda szintű hozzáférési kód szükséges, és az törli a grafikus felhasználói felületen végzett összes beállítást. Az eseménynapló nem törlődik a gyári beállítások visszaállításakor.

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

57. ábra: Fájlok és info

Ha tovább görget lefelé, megjelenik a **Készülékszoftver** gomb, ahogyan az 57. ábra jobb oldalán látható. A gomb megérintésével új készülékszoftvert telepíthet a rázógépre. Az új készülékszoftver telepítésével kapcsolatos részletes utasításokat lásd a "A készülékszoftver telepítése" a 130. oldalon.

MEGJEGYZÉS Új készülékszoftver telepítését az arra felhatalmazott szerviztechnikusnak kell elvégeznie.

<u>Szerviz</u>

A szerviz beállításokhoz való hozzáférés korlátozott és csak arra feljogosított szerviz technikusok számára megengedett.

3. 2. 4. Programok

Érintse meg a navigációs sávon a **Programok** szimbólumot a programok listájának megjelenítéséhez vagy új program létrehozásához. Lehetősége van program létrehozására, szerkesztésére, törlésére, importálására és exportálására. Az 58. ábran a programok képernyője látható új készülék (balra) és egy már egy ideje használatban lévő készülék (jobbra) esetén. A használatban lévő készülék megjeleníti a felhasználó által létrehozott programok listáját.



58. ábra: Programok

A programlistában minden bejegyzéshez tartozik egy Szem ikon (), amely megnyitja a program gyorsnézetét a program részleteivel együtt, és egy Ceruza ikon / , amely szerkesztésre nyitja meg a programot.

Előzetes beállítások létrehozása az összes programhoz

A Programok képernyőről a **Bőv. beállítások** gombbal a **Bővített beállítások** képernyőre léphet. Ezen a képernyőn határozhatja meg, hogy mi történjen egy program indításakor. Ezek a beállítások az összes létrehozott programra vonatkoznak.



59. ábra: Programbeállítások létrehozása

- 1. Érintse meg a **Bőv. beállítások** gombot.
- Válassza ki az Időzítés-vezérlés megjelenítése beállítást annak meghatározásához, hogy miként jelenjen meg a program időzítés-vezérlése:
 - » Eltelt idő: ennyi ideje fut a rázógép a Start gomb megnyomása óta, vagy
 - » Hátralevő idő: ennyi ideig fut még a rázógép, amíg le nem jár az időzítő
- Ha azt szeretné, hogy a rázógép a program indításakor azonnal elinduljon, aktiválja a Rázógép indítása a program betöltésekor beállítást. Alapértelmezés szerint program indításához a Start gombot kell megérintenie.

MEGJEGYZÉS A **Rázógép indítása a program betöltésekor** beállítás csak akkor működik az elvárásoknak megfelelően, ha zárva van a rázógép burkolata. A burkolat zárásakor a futás automatikusan elindul.

4. Nyomja meg a **Mentés** gombot.

Program létrehozása

Legfeljebb 99 programot hozhat létre és menthet el.

- 1. Nyomja meg az Új program létrehozása gombot.
- 2. Adja meg a program nevét.

Name		an to onto			_
rearrie.	10	ip to ente			_
Step 1					
Temp		Speed		Hours	Minutes
37.0	°C	50	rpm	0	1
To au	the unit	at ambient ter	mnerature k	eave the tem	n field hlank

60. ábra: Program létrehozása képernyő

 Adja meg a hőmérsékletet (kizárólag hőmérséklet-szabályozott rázógépek esetén), a fordulatszámot és a program futási idejét órákban és percekben.

Hőmérséklet-szabályozott rázógépnél a hőmérsékletszabályozás kikapcsolásához hagyja üresen a **Hőmérséklet** mezőt.

- Ha további lépést szeretne hozzáadni a programhoz, görgessen lefelé, és érintse meg a Lépés hozzáadása gombot.
- 5. Nyomja meg a Mentés gombot, a program mentéséhez.

Egy program szerkesztése

- 1. Érintse meg a ceruza ikont 🖍 a szerkeszteni kívánt program mellett.
- Módosítsa a kívánt mezőket. Nyomja meg a Mentés gombot. A programot az elvégzett módosításokkal fogja elmenteni.
- Ha lépést akar hozzáadni, a képernyő alsó szélén válassza a Lépés hozzáadása gombot.
 Ha 3 lépésnél több van, görgesse le a képernyőt ahhoz, hogy láthassa a Lépés hozzáadása gombot.

MEGJEGYZÉS Ha a több lépéses rázási folyamat közben riasztás vagy hibaüzenet jelenik meg, a rázó automatikusan leáll. Ha figyelmeztető jelzés jelenik meg, a rázási folyamat folytatódik.



61. ábra: Lépés hozzáadása

Egy program törlése

- 1. Érintse meg a ceruza ikont a szerkeszteni kívánt program mellett.
- 2. Nyomja meg a Törlés gombot. A megjelenő ablak megerősítést kér.



62. ábra: Program törlése

3. Nyomja meg a Törlés gombot. A kiválasztott program törlődik.

Egy program végrehajtása

- 1. A meglevő program végrehajtásához válassza ki a végrehajtandó programot.
- Ha a program gyorsnézetét szeretné megtekinteni, érintse meg a program melletti szem ikont
 O.
- 3. Megjeleni a Program gyors nézet ablak.



63. ábra: Program gyors nézet

4. Nyomja meg a Futtatás gombot, hogy a programot a fő képernyőre töltse.

A kiválasztott **Automatikus futás** beállításoktól függően a program azonnal, illetve a kezdőképernyő **Start** gombjának megérintése után indul el.

5. A program teljes végrehajtása után az alábbi ablak jelenik meg. Nyomja meg az OK gombot.



64. ábra: Program lezárva

Programok importálása

Módja van a rázóra olyan programot importálni, amit másik készüléken hoztak létre. Ehhez hozzá kell csatlakoztatnia az USB-meghajtót a rázógép USB-csatlakozójához.

A következő képernyő az USB-meghajtón azonosított összes programot megjeleníti.



65. ábra: Válassza ki az importálandó programot

Válassza ki azokat a programokat, amiket importálni kíván. Nyomja meg az Import gombot.

A programok sikeres importálása után az alábbi ablak jelenik meg. Nyomja meg az **OK** gombot. Az USB pendrive eltávolítható.



66. ábra: Az importálás lezárva

Programok exportálása

Lehetősége van program egyik rázóról a másikra történő exportálására. Győződjön meg arról, hogy csatlakoztattak USB pendriveot.

Ha az USB pendriveot csatolta, az alábbi képernyő jelenik meg:



67. ábra: Válassza ki az exportálandó programokat

Válassza ki azokat a programokat, amiket exportálni kíván. Nyomja meg az Export gombot.

Az exportálás lezárásakor az alábbi ablak jelenik meg. Nyomja meg az **OK** gombot. Az USB pendrive eltávolítható.



68. ábra: Exportálás lezárva

Eseménynapló

A navigációs sáv harmadik ikonja az eseménynapló, ez tartalmazza a felhasználókkal és a rendszerrel kapcsolatos eseményekről szóló adatsort Az Eseménynapló képernyő akkor jelenik meg, ha a navigációs sávon megérinti az Eseménynapló ikont a 69. ábran látható módon.



69. ábra: Eseménynapló

Ez a képernyő az aktuális események listáját jeleníti meg, dátummal és időponttal együtt.

Az egyes eseményekre vonatkozó további információk a listában található elem kiválasztásával jeleníthetők meg.

Az események a következő kategóriák szerint szűrhetők:

- Riasztások
- Figyelmeztetések
- A beállítások módosítása
- Elindítási és leállítási folyamatok
- Programfutások
- A burkolat nyílásai (kizárólag hőmérséklet-szabályozott rázógépek esetén)

Szűrő kiválasztásakor a jobb oldali Összes megjelenítése gomb BEKAPCSOLT szűrőre változik a 70. ábran látható módon.



70. ábra: Szűrt eseménynapló

Eseménynapló exportálása

 Válassza ki az exportálandó eseményeket az Exportálandó események legördülő menüből. Válassza ki a napló vagy a jelentés exportálásának formátumát.



71. ábra: Exportálandó eseménynapló kiválasztása

2. Választható egy előre beállított vagy ügyfélspecifikus dátumtartomány.

MEGJEGYZÉS Csak az elmúlt hat hónap eseményeit exportálhatja.

 Az exportálás USB pendrivera történik. A napló vagy jelentés mentéséhez USB pendriveot kell csatolni. A napló vagy jelentés letöltéséhez érintse meg a Napló exportálása gombot.



72. ábra: Csatoljon USB pendriveot az exportáláshoz

<u>Grafikonok</u>

A diagramok a fordulatszám vagy a hőmérséklet időbeli adatait (kizárólag hőmérséklet-szabályozott rázógépek esetén) mutatják. Az X tengely az időt, az Y tengely pedig a fordulatszámot vagy a hőmérsékletet mutatja.

A hőmérsékleti diagram lehetővé teszi a névleges, a tényleges és a környezeti hőmérséklet időbeli rögzítését. A **Szerkesztés** gombbal kiválaszthatja, hogy a három görbe közül melyik szerepeljen a hőmérsékleti diagramon. Ezenkívül a megjelenítési tartomány módosítható 1 napra, 7 napra vagy a napok felhasználó által meghatározott sorozatára, akár 24 teljes órára, vagy az egyes napok kiválasztott óráira.

A **Frissítés** gomb a beállításainak megfelelően frissíti a diagram megjelenítését, és megjeleníti a diagram legutóbbi betöltése vagy frissítése óta rögzített új adatokat.

Showing: 5	5/12/15, BAM to	4PM	-	3	stpoint	Actual	*	Showing: 5/1	2/15, 8AM to -	PM	Set	point	Actual	Ambie
350rpm							*	70.0°C						
300rpm							¥0	60.0°C						
250rpm								50.0°C						
200rpm								40.0°C						
150rpm								30.0°C		-			1 1	-
100rpm							Ē	20.0°C					-	
50rpm							1000	10.0°C						
Orpm	8AM 5/12/15	9AM	10AM	11AM	12PM			0.0°C L	8AM 5/12/15	9AM	10AM	11AM	12PM	

73. ábra: Fordulatszám- és hőmérsékleti diagramok

A fordulatszám- vagy hőmérsékleti diagram adatainak letöltéséhez hozzácsatlakoztathat egy USBmeghajtót a rázógép USB-csatlakozójához, és a **Diagramadatok exportálása** elem megérintésével letöltheti a fordulatszám- vagy hőmérsékleti diagram adatait.

Diagram szerkesztése

1. Érintse meg a Szerkesztés gombot a diagram szerkesztésének megkezdéséhez.



74. ábra: Diagram szerkesztése: Válassza ki a hőmérsékletet vagy a fordulatszámot a rögzítéshez

- Érintse meg a Hőmérséklet vagy a Fordulatszám gombot a megjelenítendő diagram kiválasztásához.
- 3. Görgessen tovább lefelé a Dátumtartomány beállításainak megjelenítéséhez.
- Érintse meg az 1 nap, 7 nap vagy a Felhasználó által meghatározott lehetőséget egy egyéni dátumtartomány kiválasztásához.
- Forgassa el a beállítókerék forgókapcsolóját a kezdődátum kiválasztásához, és érintse meg az Ettől gombot.

 Forgassa el a beállítókerék forgókapcsolóját a záródátum kiválasztásához, és érintse meg az Eddig gombot.



75. ábra: Dátumtartomány kiválasztása a diagramhoz

- 7. Görgessen tovább lefelé az Időtartomány beállítások megjelenítéséhez.
- Érintse meg a napi 24 óra gombot, vagy érintse meg a Felhasználó által meghatározott gombot egy adott időszak megjelenítéséhez, és
 - Forgassa el a beállítókerék forgókapcsolóját a napi kezdési időpont kiválasztásához, és érintse meg a Kezdet gombot.
 - Forgassa el a beállítókerék forgókapcsolóját a napi befejezési idő kiválasztásához, és érintse meg a **Befejezés** gombot.

MEGJEGYZÉS A kezdő és befejező időpontok kiválasztásával dönthet a két időpont közötti időszak pontos ábrázolásáról. Ha a napi 24 órás lehetőséggel dolgozik, a rögzítés minden nap 00:00 órakor kezdődik és 23:59 órakor fejeződik be.



76. ábra: Napszakok kiválasztása a rögzítéshez

9. A módosítások vagy az ügyfélspecifikus beállítások mentéséhez érintse meg a Mentés gombot.

3.3. Tartozékok



Sérülésveszély és biológiai veszélyeztetés a törött edények miatt.

A szakszerűtlenül felszerelt tartozékok miatt az üvegek eltörhetnek és a minták kiömölhetnek.

VIGYÁZAT

Biztosítsa, hogy a tartozékok megfelelően legyenek felszerelve, a megfelelő szerszámok és csavarok használatával.

Biztosítsa, hogy a tartozékok megfelelően rögzüljenek a tálcára. Az edényeket mindig csak az adott méretnek megfelelő tartozékokkal együtt használja.



VIGYÁZAT

Az éles szélek vágott sérüléseket okozhatnak.

Óvatosan bánjon a tálcával és más tartozékokkal.



A mozgó tálca becsípheti az ujjait.

Sose tegye az ujjait a mozgó tálcára vagy az alá.

VIGYÁZAT

MEGJEGYZÉS Az üzemeltető felelős a tartozékok szabályos felszereléséért.

A telepítés itt leírt eljárása további intézkedésig a rázó kezelési útmutatóban felsorolt összes kivitelére érvényes.

Mindig a megfelelő szerszámokat használja, különösen a tartozékokhoz mellékelteket. Ha ezek közül néhány elveszett, újrarendelheti a felsorolt pótalkatrészkészleteket ("Tartozékok" a 17. oldalon). Kizárólag a megadott szerszámot és csak a mellékelt csavarokat használja.

3. 3. 1. A tálca felszerelése

VIGYÁZAT Az egyes rázógépek kompatibilis tálcáinak teljes listáját lásd a "Tálcák" a 18. oldalon. A hőmérséklet-szabályozott rázógépek kizárólag a gyárilag előre beszerelt tálcákkal üzemeltethetők.

VIGYÁZAT Az éles szélek vágott sérüléseket okozhatnak. Tálca kiszerelésekor nyúljon a tálca alá. Sose emelje meg a tálcát a rászerelt lombiktartónál fogva. Óvatosan bánjon a tálcával és más tartozékokkal.

VIGYÁZAT A Solaris 4000 I/4000 R rázógép 11 x 14-es tálcájának mozgása közben fennáll az ujjak és a kéz becsípődésének a veszélye. Ne használja a 11 x 14-es tálcát Solaris 4000 I/4000 R rázógépekhez!

VIGYÁZAT Kizárólag a tálcával szállított rögzítő szerszámokat és csavarokat használja. Más csavarok vagy nem megfelelő rögzítő szerszámok használata nem megfelelő telepítést jelent és ezzel a rázó és tartozékainak esetleges károsodását okozza.

A tálcát mindig az összes csavarjával építse be.

Mindig a megfelelő rögzítő szerszámot használja:

- 3/16 colos rögzítőszerszám (GT530066) a Solaris 2000, Solaris 2000 I/R és 4000 I/R rázógépek összes tálcájához
- 7/32 colos rögzítőszerszám (GT530080) a Solaris 4000 rázógépek összes tálcájához

Univerzális tálcák

Valamennyi Solaris 2000 és 4000 sorozatú rázógépet univerzális tálcával, csavarokkal és rögzítőszerszámmal szállítunk. Alkalmazásaihoz további tálcákat vásárolhat. Az egyes rázógépek kompatibilis tálcáinak teljes listáját lásd az "Tálcák" a 18. oldalon.

A 2000 I, 2000 R, 4000 I és 4000 R rázógépeknél a tálca gyárilag be van építve a rázógépbe. Sem a tálca, sem a csavarok nem kaphatók külön alkatrészként.



78. ábra: Univerzális tálca egy Solaris 2000 (bal) és egy Solaris 4000 (jobb) rázóra szerelve

1. Óvatosan helyezze a tálcát vízszintesen a rázó 3 szerelési pontjára.

A Solaris 2000 rázógépek tálcái 3 rögzítőfurattal rendelkeznek. A Solaris 4000 rázógépek tálcái 4 rögzítőfurattal rendelkeznek.

A Solaris 2000 I/R és 4000 I/R rázógépek tálcái 3 rögzítőfurattal rendelkeznek.

 A tálca rázóra való rögzítéséhez óvatosan húzza meg a csavarokat. Ha a rögzítő szerszám kezd meggörbülni, ne húzza tovább a csavart.

Kétlépcsős tálcák

MEGJEGYZÉS A kétlépcsős tálcák nem alkalmasak Solaris 2000 I/R és 4000 I/R rázógépekhez.



79. ábra: Kétlépcsős tálcaegység

Kapcsolja össze a felső és alsó tálcát, a 4 sarkába szerelt támasztókkal. A támasztókat a megfelelő csavarokkal rögzítse a felső és az alsó tálcához.

A támasztó tálcához való rögzítéséhez óvatosan húzza meg a csavarokat. Ha a rögzítő szerszám kezd meggörbülni, ne húzza tovább a csavart.

Az alsó tálcán vannak a rázóhoz történő rögzítést szolgáló szerelő furatok. Biztosítsa, hogy az összeszereléskor az alsó tálca legyen az alsó oldalon.

Tartozékok

MEGJEGYZÉS Mielőtt az edényeket a tálcára helyezi, még egyszer ellenőrizze, hogy nem lötyög-e a tálca szerelvény.



80. ábra: Kétlépcsős tálca a Solaris 4000 rázón

1. Óvatosan helyezze a kétlépcsős tálcát vízszintesen a rázó szerelő lapjára.

A Solaris 2000 tálcáin 3, a Solaris 4000 tálcáin 4 szerelési furat található.

 A tálca rázóra való rögzítéséhez óvatosan húzza meg a csavarokat. Ha a rögzítő szerszám kezd meggörbülni, ne húzza tovább a csavart.

3. 3. 2. A dugattyútartó és az edény behelyezése

VIGYÁZAT A törött vagy nem záró edények biológiai veszélyt jelentenek. A szakszerűtlenül felszerelt tartozékok miatt a minták kiömölhetnek. Biztosítsa, hogy a tartozékok megfelelően legyenek felszerelve, a megfelelő szerszámok és csavarok használatával. A behelyezéshez 150 mm/6 col szárhosszal rendelkező csavarhúzó (PH2) szükséges (cikkszám: 75004131). Biztosítsa, hogy a tartozékok megfelelően rögzüljenek a tálcára. Az edényeket mindig a méretükhöz illő tartozékokkal használja. Az edények legyenek épek és megfelelően felszerelve.

VIGYÁZAT Az éles szélek vágott sérüléseket okozhatnak. Óvatosan bánjon a tálcával és más tartozékokkal.

VIGYÁZAT A dugattyútartó biztonságos rögzítése csak a mellékelt csavarokkal együtt garantált. Helyezze be a mellékelt csavarokat a lombiktartóban található összes rögzítőfuratba.

MEGJEGYZÉS Ha a csavarok túl hosszúak, csökkenhet a cserélhető tálca stabilitása és funkcionalitása. Ügyeljen arra, hogy csak a mellékelt csavarokat használja a rögzítéshez.

Lombiktartó

Minden lombiktartó egy edénytartókapocsból, a lombik/palack nagyságától függően egy vagy két tartórugóból és csavarokból áll, amik a tálcához rögzítik. Mindig a lombiktartóval szállított csavarokat használja.



① Rugós hüvely ② Rugó ③ Csavarok ④ Tartókapocs ⑤ Rugó

81. ábra: Lombiktartó 2 rugóval

- 1. Ha szükséges, rögzítse a rugót a tartókapcsokhoz az ábrákon látható módon.
- A rugós hüvelyeket az ábrán látható módon kell a tartókapcsok közé tolni. Egyes lombiktartók 2 rugót tartalmaznak. A második rugót a tartókapocs alsó végéhez szerelik, miután rögzítették a lombiktartót a tálcára.
- 3. A vele szállított csavarokkal rögzítse a tartókapcsokat a tálcára.

Edény

- Óvatosan helyezze be a kívánt edényt a lombiktartóba. Ehhez előbb húzza el a rögzítőrugót annyira, hogy az edény alja a lombiktartó közepére kerüljön. Hagyja az edényt lazán a helyre siklani és rögzítse a lombiktartó szélesebb fenekénél. A rugó biztos helyzetben tartja az edénynyakot.
- Mielőtt a rázót használatba veszi, győződjön meg arról, hogy az összes edényt megfelelően rögzítették.

Lehetőség szerint az edényt dugóval kell lezárni, hogy a keverési folyamat közben ne kerüljenek ki anyagok.

3. 3. 3. Szögletes lombiktartó felszerelése

VIGYÁZAT A törött vagy nem záró edények biológiai veszélyt jelentenek. A szakszerűtlenül felszerelt tartozékok miatt a minták kiömölhetnek. Biztosítsa, hogy a tartozékok megfelelően legyenek felszerelve, a megfelelő szerszámok és csavarok használatával. A behelyezéshez 150 mm/6 col szárhosszal rendelkező csavarhúzó (PH2) szükséges (cikkszám: 75004131). Biztosítsa, hogy a tartozékok megfelelően rögzüljenek a tálcára. Az edényeket mindig a méretükhöz illő tartozékokkal használja. Az edények legyenek épek és megfelelően felszerelve.

VIGYÁZAT Az éles szélek vágott sérüléseket okozhatnak. Óvatosan bánjon a tálcával és más tartozékokkal.



Tartozékok

Szögletes lombiktartó

Valamennyi szögletes lombiktartó egy rögzítőfuratokkal rendelkező alaplemezből és négy rugós tartókapocsból áll. Tartalmazza továbbá a csavarokat, amelyek a lombiktartó tálcához történő rögzítéséhez szükségesek. Mindig a lombiktartóval szállított csavarokat használja.



① Rögzítőfuratok ② Rugó ③ Tartókapocs ④ Csavarok

82. ábra: Szögletes lombiktartó

- 1. Szükség esetén akassza be az összes rugót 2 a tartókapcsokba 3 a 82. ábran látható módon.
- 2. Helyezze a tálcára a rögzítőegységet, és igazítsa be a rögzítőfuratokat ①.
- 3. Erősítse hozzá a rögzítőegységet a tálcához a mellékelt csavarokkal ④.
- 4. Helyezze be felülről a négyzet alakú edényt.
- Győződjön meg arról, hogy az edény felfekszik az alaplapra, és stabilan helyezkedik el a lombiktartóban.

3. 3. 4. A kémcsőállvány felszerelése

VIGYÁZAT Az éles szélek vágott sérüléseket okozhatnak. Óvatosan bánjon a tálcával és más tartozékokkal.

Minden kémcsőállvány két rögzítőkengyelből (drótkeret vagy lemez) áll, amelyekhez egy-egy rögzítőcsap-fej és kémcsőállványonként két recés fejű csavar tartozik, a tálcához történő rögzítés céljából. Minden kémcsőállvány részegységet előre összeszerelve szállítunk, kémcsőállvánnyal és habbetéttel (habbetétekkel) együtt. Mindig az állványhoz mellékelt csavarokat használja.



83. ábra: Kémcsőállvány részegység rögzítőkonzolokkal (drótkeret)



Kémcsőállvány
 Szivacs betétek
 Rögzítőcsapos fej
 Szerelőkonzol

84. ábra: Kémcsőállvány részegység rögzítőkonzolokkal (lemez)

- Szerelje fel a szerelőkonzolokat a tálcára a recés fejű csavarokkal. A recézett csavarokat kézzel húzza meg.
- 2. Helyezze be a kémcsőállványt a rögzítőkonzolokba.
- A rögzítőcsapos fejekkel állítsa be a megfelelő szöget. Kézzel húzza meg a rögzítőcsapos fejeket.

3. 3. 5. Állítható szögű kémcsőállvány-tartó felszerelése

VIGYÁZAT Az éles szélek vágott sérüléseket okozhatnak. Óvatosan bánjon a tálcával és más tartozékokkal. A kémcsőállvány-tartók és kémcsőállványok különböző méretekben is elérhetők. Egy felsorolás található az "Állítható szögű kémcsőállványok" a 28. oldalon. Valamennyi kémcsőállvány-tartó hét pozícióban állítható, mindkét irányban 0°-tól elforgatható, és 15°, 30° és 45°-ban rögzíthető. Mindig az állványtartóhoz mellékelt csavarokat használja.



85. ábra: Állítható szögű kémcsőállvány-tartók kémcsőállvánnyal felszerelve

- 1. Szerelje fel az állványtartót ① a rázógép tálcájára a mellékelt csavarok segítségével.
- Feszítse szét finoman a fémfüleket ③ az állványtartó forgatható részének ⑥ mindkét végén, és helyezze be a kémcsőállványt ②.
- 3. Rögzítse a kémcsőállványt 2 a fülek óvatos meglazításával 3.
- Húzza ki a rögzítőcsapok fejét (5) a tartó mindkét végén. A csapok a fej 1/4 fordulatnyi elfordításával reteszelhetők ki.
- 5. Fordítsa el az állványtartó 6 forgatható részét 1 a kívánt 15°, 30° vagy 45°-os szögbe.
- Helyezze vissza a rögzítőcsapot (5) a pozicionáló furatba (4), és rögzítse a csapot a fej 1/4 fordulattal történő elfordításával.
- Az állvány eltávolításához feszítse szét a fémfüleket ③ a forgatható rész mindkét végén ⑥, és emelje ki a kémcsőállványt ①.

3. 3. 6. Lombiktartó felszerelése Mikrotiter/Deepwell tálcákhoz

VIGYÁZAT Az éles szélek vágott sérüléseket okozhatnak. Óvatosan bánjon a tálcával és más tartozékokkal.



- Deepwell-tálca
- 2 Csavar
- ③ A Mikrotiter-tálca lombiktartója

86. ábra: Lombiktartó részegység Mikrotiter/Deepwell tálcákhoz

- 1. Helyezze a Mikrotiter-tálca keretét a tálcára.
- A Mikrotiter- / Deepwell-tálcakészlettel szállított csavarok segítségével szerelje a Mikrotitertálca keretét a tálcára.
- 3. Illessze a Mikrotiter- vagy Deepwell-tálcát a Mikrotiter-tálca keretébe.
- 4. Az óvatos megemelésével győződjön meg a Mikrotiter- vagy Deepwell-tálca szilárd rögzítéséről.

3. 3. 7. Állítható edénytartó felszerelése

VIGYÁZAT Az éles szélek vágott sérüléseket okozhatnak. Óvatosan bánjon a tálcával és más tartozékokkal.





[®] Csavarok

87. ábra: Egy fokozatban állítható magasságú edénytartó részegység

Tartozékok





① Rögzítőfuratok ② Állítható rudak ③ Rögzítőcsap-fejek ④ A tároló oldalfala ⑤ Csúszásgátló szőnyeg
 ⑥ Csavarok

88. ábra: Két fokozatban állítható magasságú edénytartó részegység

- Emelje fel a csúszásgátló szőnyeget, hogy láthatóvá váljanak a rögzítőfuratok (tétel ① a 88. ábran).
- 2. Helyezze a tálcára az állítható edénytartót, és igazítsa be a rögzítőfuratokat.
- 3. Rögzítse hozzá a tálcához az állítható edénytartót a mellékelt csavarokkal 6.
- 4. Helyezze a csúszásgátló szőnyeget 5 az állítható edénytartó aljára.
- 5. Helyezze az edény(eke)t a csúszásgátló szőnyegre ⑤.
- 6. Lazítsa meg a rögzítőcsapok fejét 3.
- 7. Állítsa be az állítható rudak @ helyzetét úgy, hogy az edényeket biztonságosan rögzüljenek.
- 8. Húzza meg újra a rögzítőcsapok 3 fejét.
- 9. Óvatosan felemelve győződjön meg arról, hogy az edény biztonságosan van rögzítve.

3. 3. 8. A főzőpohár állványok felszerelése

VIGYÁZAT Az éles szélek vágott sérüléseket okozhatnak. Óvatosan bánjon a tálcával és más tartozékokkal.



89. ábra: Főzőpohár állvány részegység

- 1. Helyezze az alsó lemezt 2 a tálcára.
- 2. Helyezze a főzőpohár állványt ① az alsó lemezre.
- 3. Igazítsa hozzá a tálcához a rögzítőfuratokat.
- 4. Rögzítse a főzőpohár állvány részegységet a mellékelt csavarokkal ③.

3. 3. 9. A szétválasztó tölcsértartó felszerelése

VIGYÁZAT Az éles szélek vágott sérüléseket okozhatnak. Óvatosan bánjon a tálcával és más tartozékokkal.

MEGJEGYZÉS Ügyeljen arra, hogy a megfelelő csavarokat használja. A csomag kétféle csavart tartalmaz. A hosszabb, piros menetzárral ellátott csavarokat a függőleges állvány rögzítéséhez kell használni. A rövidebb, menetzár nélküli csavarok kizárólag a szétválasztó tölcsértartó tálcára szereléséhez használhatók.

A szétválasztó tölcsértartó alapvetően kétféle helyzetben szerelhető fel a rázógép tálcájára: vízszintesen és függőlegesen.



90. ábra: A szétválasztó tölcsértartó beszerelési helyei

A 90. ábran látható ③ vízszintes helyzet olyan magasságban állítható be, hogy a ④ szerinti ferde beigazítást kapjuk. Ezen túlmenően a szétválasztó tölcsértartó a (tartozékként kapható) függőleges állvánnyal is kombinálható, így létrehozhatók az ① és ② ábrákon látható függőleges elrendezések. A tartó felszerelhető úgy is, hogy a tölcsércsap lefelé nézzen ①, vagy fordítva, így a ③ ábrán látható elrendezést kapjuk.

Függőleges beszerelés előkészítése



91. ábra: Szétválasztó tölcsértartó (balra) függőleges állványhoz való tartozékokkal (jobbra)

- Rögzítse hozzá a szétválasztó tölcsértartót a függőleges állvány hosszú végéhez az ① és
 ② szerint, a 90. ábran látható módon.
- Igazítsa hozzá azt a függőleges állvány tölcsértartójához tartozó négy rögzítőfurathoz (@ a 91. ábran) az alábbiak szerint:
 - A tölcséres csapot a szabványos függőleges beszereléshez használt furatokkal (③ a 91. ábran) lefelé fordítva kell beszerelni.
 - Alkalmazza a függőlegesen elfordított beszereléshez használt furatokat (@ a 91. ábran) a tölcséres csap felfelé néző beszereléséhez.
 - c. Rögzítse hozzá a szétválasztó tölcsértartót a függőleges állványhoz a mellékelt csavarokkal (⑤ a 91. ábran).

Felszerelés a tálcára

- 1. Helyezze a szétválasztó tölcsértartó részegységet a tálcára.
- 2. Igazítsa hozzá a tálcához a rögzítőfuratokat.

A tölcsér felszerelése



① Szétválasztó tölcsér
② Hátsó tölcsértartó
③ Elülső tölcsértartó
④ Rögzítőcsap fej, magasságállítás
⑤ Rögzítő csapfej, hosszbeállítás
⑥ Menetzár nélküli csavarok

92. ábra: Szétválasztó tölcsértartó részegység

- Lazítsa meg a hosszbeállítás rögzítő csapfejét (5), és csúsztassa el az elülső tölcsértartót (3) teljes mértékben előre.
- 2. Helyezze be a tölcsér hátsó végét ① a hátsó tölcsértartóba ②.
- Csúsztassa el hátrafelé az elülső tölcsértartót ③, és helyezze be a tölcsér nyakát az elülső tölcsértartó nyílásába ③.
- 4. Állítsa az elülső tölcsértartót ③ megfelelő magasságba, és rögzítse a magasságállító rögzítő csapfej ④ meghúzásával. A magasságállítás megváltoztatásával ferde helyzetbe állíthatja a szétválasztó tölcsért.
- 5. Rögzítse az elülső tölcsértartót ③ a hosszbeállító rögzítő csapfej ⑤ meghúzásával.

3. 3. 10. Többfunkciós tároló felszerelése

VIGYÁZAT Az éles szélek vágott sérüléseket okozhatnak. Óvatosan bánjon a tálcával és más tartozékokkal.



93. ábra: Többfunkciós tároló részegység

A telepítéshez az alábbiak szerint járjon el:

- 1. Helyezze a tálcára a többfunkciós tárolót 2.
- 2. Igazítsa hozzá a tálcához a rögzítőfuratokat.
- 3. Rögzítse a többcélú tárolót 2 a hozzá mellékelt csavarokkal 3.
- 4. Helyezze a gumiszőnyeget ① a tárolóba②.
- 5. Helyezze a palackokat a tárolóra, és rögzítse őket a mellékelt O-gyűrűkkel ④, lásd 94. ábra.

MEGJEGYZÉS 100 fordulat/percnél nagyobb fordulatszám esetén O-gyűrűk használata javasolt.



94. ábra: Többfunkciós tároló rögzített tárolókkal és O-gyűrűkkel

MEGJEGYZÉS A tartályok rögzítésekor ügyeljen arra, hogy ne feszítse meg túlzottan az O-gyűrűket. Ha minden egyes O-gyűrűt megnyújt az összeszereléshez, érez majd egy holtpontot a rugalmasság ellenőrzésekor, amelyen túl az O-gyűrűket nem lehet megnyújtani. Ha elérte ezt a pontot, ne próbálja tovább feszíteni az O-gyűrűket, mert megrepednek.

3. 3. 11. Az öntapadós szőnyeg lerakása és használata

Az öntapadós szőnyegek széles, lapos fenekű tartályokhoz, például Erlenmeyer-lombikokhoz, Fernbach-lombikokhoz, médialombikokhoz, Mikrotiter tálcákhoz, Petri-csészékhez, sejttenyésztő lombikokhoz, térfogatmérő lombikokhoz és főzőpoharakhoz alkalmasak. Általában bármilyen olyan edény használható, amely képes függőlegesen állni, de ahogy az alapterület és a magasság aránya csökken, úgy csökken a maximálisan elérhető rázási fordulatszám is. A magas, kis fenekű tartályok kevésbé alkalmasak, lásd a 95. ábrat. A nagy Erlenmeyer-lombikok (például a 3000 ml-esek) erősebben tapadnak, mint a kicsik (például a 100 ml-esek).



95. ábra: Öntapadós szőnyeg: Megfelelő lombikformák

Általában kétféle öntapadós szőnyeg különböztethető meg:

- Alacsony tapadású szőnyeg (átlátszó) 75004111 és 75004117
 - » egyszerű kezelhetőség az alacsonyabb tapadásnak köszönhetően
 - » különösen főzőpoharakhoz alkalmas
 - » bármilyen edényhez alkalmas 200 fordulat/perc maximális fordulatszámig (sejtkultúrákhoz történő alkalmazás)
- Nagy tapadású szőnyeg (fekete, átlátszó) 75004126 és 75004127
 - » megbízható edényrögzítés a nagyobb tapadásnak köszönhetően
 - » bármilyen edényhez alkalmas
 - » maximális fordulatszám 400 fordulat/perc (akár 500 ml-es Erlenmeyer-lombikhoz is)

<u>Eszközök</u>

A 96. ábran látható felhordó tekercs az öntapadós szőnyegek tartozékaként kerül forgalomba. Segédeszköz, amely megkönnyíti az öntapadós szőnyeg tálcára nyomását.



96. ábra: Felhordó tekercs öntapadós szőnyegekhez

Száraz szőnyeg behelyezése

- 1. Győződjön meg arról, hogy a rázógép tálcája tiszta és száraz.
- 2. Távolítsa el a védőfóliát az öntapadós szőnyeg egyik oldaláról.
- 3. Ragassza hozzá erősen a szőnyeget a tálcához a ragasztós oldallal lefelé.
- 4. Távolítsa el a védőfóliát a felső oldalról.

Nedves szőnyeg behelyezése

- 1. Távolítsa el a védőfóliát az öntapadós szőnyeg mindkét oldaláról.
- 2. Áztassa be vízbe a szőnyeget, vagy alaposan öblítse át.
- 3. Hagyja lecsöpögni a vizet a szőnyegről.
- 4. Terítse a nedves szőnyeget a tiszta tálcára, és helyezze a kívánt pozícióba.
- 5. Használat előtt hagyja a szőnyeget 24 órán át száradni a tálcán.

Edényjellemzőkkel kapcsolatos megjegyzések

A maximális fordulatszám az edény típusától függ, amely méret, alak, fenékforma, anyag, felületi érdesség és tisztaság szempontjából eltérő. A rázógép használata előtt ellenőrizze ezeket a paramétereket, hogy meghatározhassa az adott alkalmazáshoz ideális szőnyeget.

Tekintse meg az 53. táblázatban, hogy a laboratóriumi anyag alkalmas-e az öntapadós szőnyegekkel történő együttes használatra.

Műanyag-/üvegtermékek	A tapadás értékelése	Fém	A tapadás értékelése
Üveg (boroszilikát üveg)	5	Acél, rozsdamentes	5
Polipropilén (PP)	3	Alumínium	5
Polipropilén kopolimer (PPCO)	3	Eloxált alumínium	4
Polietilén (HPDE) (LPDE)	3		
Polikarbonát (PC)	4		
Polietilén-tereftalát (PET)	4		
Polietilén-tereftalát-kopoliészter	4		
(PETG)			
Polimetilpentén (PMP)	3		
Politetrafluor-etilén (PTFE)	0*		
Tetrafluor-etilén-hexafluor-	0*		
propilén kopolimer (FEP)			
Perfluor-alkoxi polimerek (PFA)	2		
Szilikon	1*		
A			

Értékelések: 0 = nincs tapadás; 5 = legjobb tapadás/*ne használjon 0 és 1 értékelésű edényeket (PTFE, FEP és szilikon).

53. táblázat: Az edény anyagainak kompatibilitása

MEGJEGYZÉS Kérjük, vegye figyelembe, hogy a maximális rázási fordulatszám csak az 53. táblázatban szereplő, 5-ös értékelésű anyagokkal érhető el.

MEGJEGYZÉS Az üvegtől eltérő anyagok esetében Önnek kell kipróbálnia és meghatároznia, hogy melyik anyagtípus milyen terheléshez, fordulatszámhoz, futási időhöz és hőmérséklethez alkalmas.
Fordulatszám-tartomány és futási idő korlátozás üvegedényekhez (30%-os töltöttség)

Üvegedények fordulatszám és futási idő határértékei szobahőmérsékleten (30%-os töltöttség). A magasabb hőmérséklet csökkenti a tapadást. Az edények a hőmérséklet megváltoztatásával leválhatnak a szőnyegről. Kezdje a rázási folyamatot akklimatizált edényekkel.

	Mikrotiter/Deepwell	Főzőpohár			
	96 x 2 ml	25 ml	125 ml	250 ml	500 ml
125 ford./perc	nincs korlátozva a futási id	lő			
150 ford./perc	72 óra	72 óra	72 óra	72 óra	72 óra
200 ford./perc	72 óra	72 óra	72 óra	72 óra	72 óra
250 ford./perc	72 óra				
300 ford./perc					
350 ford./perc					
400 ford./perc					

54. táblázat: Alacsony tapadású szőnyeg – a fordulatszám-tartomány és futási idő korlátozása szobahőmérsékleten

	Mikrotiter/ Deepwell	Erlenme	yer						
	D96 x 2 ml	25 ml	125 ml	250 ml	500 ml	1000 ml	2000 ml	3000 ml	5000 ml
125 ford./	nincs korlátozva	a a futási io	dő						
perc									
150 ford./	72 óra	72 óra	72 óra	72 óra	72 óra	72 óra	72 óra	72 óra	72 óra
perc									
200 ford./	72 óra	72 óra	72 óra	72 óra	72 óra	72 óra	72 óra	72 óra	72 óra
perc									
250 ford./	72 óra	72 óra	72 óra	72 óra	72 óra	72 óra	72 óra	48 óra	24 óra
perc									
300 ford./	72 óra	72 óra	72 óra	72 óra	72 óra	72 óra	48 óra	24 óra	
perc									
350 ford./	48 óra	48 óra	48 óra	48 óra	48 óra	48 óra	24 óra		
perc									
400 ford./	24 óra	24 óra	24 óra	24 óra	24 óra				
perc									

55. táblázat: Nagy tapadású szőnyeg – a fordulatszám-tartomány és futási idő korlátozása szobahőmérsékleten

<u>Előkészítés</u>

VIGYÁZAT Üvegtermékek kezelésekor mindig viseljen védőszemüveget és kesztyűt.

Mielőtt az edényeket az öntapadós szőnyegre helyezné:

- Ellenőrizze az edényeket az esetleges károsodások szempontjából. Ne használjon károsodott edényeket!
- 2. Törölje át az edényt tiszta és száraz ruhával a nedvesség teljes eltávolításához.
- 3. Távolítsa el a zsírt és egyéb szennyeződéseket.
- 4. Ellenőrizze, hogy az edény alja és az öntapadós szőnyeg száraz és tiszta-e.

VIGYÁZAT Már egy alig látható vékony réteg vagy egyetlen vízcsepp is az edény alján azt okozhatja, hogy az edény egy bizonyos futási idő után elveszíti a tapadását és leválik a szőnyegről.



97. ábra: Öntapadós szőnyeg: törölje szárazra az edényt

5. A kiömlés elkerülése érdekében csökkentse a töltési szinteket, vagy zárja le az edényeket.

VIGYÁZAT Ha az edény tökéletesen is tapad, a külső falon végigfutó egyetlen csepp folyadék is azt okozhatja, hogy az edény egy bizonyos futási idő után leválik a szőnyegről.

Edények behelyezése

- 1. Helyezze az edényeket a szőnyegre, és nyomja meg lefelé azokat.
- 2. Óvatosan nyomja vagy húzza őket, hogy az összes edény megbízhatóan tapadjon a szőnyegre.

MEGJEGYZÉS Vegye figyelembe, hogy alacsony hőmérsékleten vagy hosszan tartó működés közben kondenzvíz képződhet, ami a tartályok nem kívánt leválásához vezethet. A minta és az edény hőmérséklete nem térhet el a rázógép megadott hőmérséklettartományától. A rázógép lehetséges hőmérsékleti tartományára vonatkozó további információk a "Műszaki adatok" a 11. oldalon című részben találhatók.

Tartozékok

Edények eltávolítása

- 1. Érintse meg a kijelzőn a Stop gombot, vagy várja meg, amíg befejeződik a futás.
- 2. Várja meg, amíg a tálca teljesen megáll.
- Óvatosan és egyenletesen húzza az edény nyakát oldalra, és várjon néhány másodpercet. A nagy fenékfelülettel rendelkező edények esetében a leválás néhány másodpercig is eltarthat.



98. ábra: Öntapadós szőnyeg: Óvatosan tolja vagy húzza le az edényt

VIGYÁZAT Óvatosan és nagy erő alkalmazása nélkül nyomja vagy húzza! Az edények hirtelen meglazulhatnak, és kiömölhet a tartalmuk!

4. Ha az edényt nem lehet meglazítani, öntsön egy kis vizet az aljára.

Hagyja a szőnyeget megszáradni a levegőn, mielőtt újra használná azt.

A szőnyeg eltávolítása

 Az öntapadós szőnyeg leválasztásához emelje meg az egyik oldalát, és lassan húzza el a tálcától a 99. ábran látható módon.



99. ábra: Öntapadós szőnyeg: Emelje fel és húzza el a szőnyeget az eltávolításhoz

Tisztítás és fertőtlenítés

Ha csökken a tapadóképesség, tisztítsa meg a szőnyeget.

VIGYÁZAT Ne használjon semmilyen oldószert, mosószert vagy tisztítószert, kivéve bőrápoló adalékanyagokat nem tartalmazó, enyhe mosogatószappant. Az ajánlott koncentráció max. 2 ml tisztítószer 1 liter vízhez.

Tartozékok

MEGJEGYZÉS A szőnyeg közvetlen érintkezése a koncentrált tisztítószerrel csökkenti a tapadást. Ebben az esetben az öblítés önmagában nem elegendő. Ha közvetlenül tisztítószereket használ, öblítse le a szőnyeget, áztassa vízben 24 órán át, majd szárítsa 24 órán át a teljes tapadás helyreállítása érdekében.

MEGJEGYZÉS Kizárólag 70%-os izopropanollal vagy 70%-os etanollal fertőtlenítsen.

Enyhén szennyezett vagy poros:

- 1. A tálcán lévő szőnyegeket nedves ruhával felülről törölje át, vagy fertőtlenítse.
- 2. Nedvesítse meg a szőnyeget.
- 3. Törölje le a maradványokat szivaccsal.
- 4. Törölje át a szőnyeget tiszta vízzel átitatott szivaccsal.
- 5. Szükség esetén fertőtlenítse a szőnyeget.
- 6. Hagyja a szőnyeget 24 órán át száradni.

Erősen szennyezett:

- 1. Vegye le a szőnyeget a tálcáról, hogy teljesen megtisztítsa azt.
- 2. Áztassa be a szőnyeget hígított mosogatószerrel kevert vízbe.
- 3. Alaposan dörzsölje át a szőnyeget egy szivaccsal, hogy minden maradványt eltávolítson.
- 4. Tisztítás után alaposan öblítse le a szőnyeget vízzel.
- 5. Szükség esetén fertőtlenítse a szőnyeget.
- 6. Hagyja a szőnyeget 24 órán át száradni.
- A behelyezés a "Száraz szőnyeg behelyezése" a 107. oldalon vagy a "Nedves szőnyeg behelyezése" a 107. oldalon leírtak szerint történik.

3. 3. 12. Gázelosztó felszerelése

A gázelosztó lehetővé teszi inert gáz, például nitrogén vagy szén-dioxid bevezetését több edénybe a rázógép működése közben (rendelési információkért lásd az "Általános tartozékok" a 42. oldalon).

A gázelosztó a Solaris 2000 I/R és 4000 I/R rázógépek gyárilag beszerelt felszerelése, azonban vevő által beszerelhető tartozékként is elérhető, ha eredetileg nem volt felszerelve a készülékre. Valamennyi Solaris 2000 I/R és 4000 I/R rázógépnél a két oldalsó panel egyikén található egy tömszelence (a helyére vonatkozó információkat lásd "Termék áttekintés" a 44. oldalon) a laboratóriumi gázellátó tömlő elosztóegységhez való vezetéséhez. A gázelosztónak egy bemeneti nyílása van a gázellátáshoz, és nyolc kimeneti nyílása a tartályok számára.



- 1 Laboratóriumi gázellátó tömlő
- ② Gázelosztó részegység
- ③ Tömlős kivezetések
- ④ Beömlőnyílás
- 5 Rögzítőcsavarok

100. ábra: Solaris 4000 R rázógépbe szerelt gázelosztó

A telepítéshez az alábbiak szerint járjon el:

- 1. Rögzítse hozzá a gázelosztót 2 a rázókamra hátsó falához a mellékelt csavarokkal 5.
- Távolítsa el a műanyag dugót a tömszelence tömítésének külső oldaláról (a helyére vonatkozó információkat lásd az "Termék áttekintés" a 44. oldalon).
- 3. Vágókéssel vágjon keresztet a rázókamra belsejében található gumitömlő membránjába.
- 4. Vágjon le egy tömlődarabot a laboratóriumi gázellátó tömlőnek megfelelő hosszúságúra ①.
- Csatlakoztassa hozzá a laboratóriumi gázellátó tömlőt ① az elosztó bemeneti nyílásának kampójához ④.
- 6. Győződjön meg arról, hogy a laboratóriumi gázellátó tömlő ① nem ér hozzá a burkolat karjaihoz, és vezesse ki a tömszelence gumitömítésén keresztül.

Tartozékok

- 7. Csatlakoztassa hozzá a laboratóriumi gázellátó tömlőt ① a gázellátáshoz.
- Vágjon le 8 db tömlődarabot, és csatlakoztassa hozzá őket az elosztó ③ kivezetéseihez. 8-nál több edény esetén használjon Y-csatlakozókat; 8-nál kevesebb edény esetén a nem használt tömlőket válassza le a gázveszteség csökkentése érdekében.
- 9. Használjon megfelelő dugókat és tömítéseket a csövek edényekbe vezetéséhez.

3. 4. Feltöltés és rendeltetésszerű használat



FIGYEL-

MEZTETÉS

A kiváltott kémiai reakciók tüzet okozhatnak.

Ne kapcsolja be a rázót olyan fordulatszámmal, aminél az edények tartalma kiömölhet.

Csak lassan emelje a fordulatszámot. Vegyszerekkel történő használat előtt javasoljuk, hogy végezzen próbát vízzel.



A szakszerűtlen feltöltés és a sérült tartozékok befolyásolhatják a biztonságot.

VIGYÁZAT

- Biztosítsa, hogy a töltet (minták és tartozékok) a tálca közepéhez képest szimmetrikusan legyen elosztva. Kétlépcsős tálca használata esetén ügyeljen arra, hogy az alsó tálca jobban meg legyen rakva, mint a felső.
- Soha ne terhelje túl a rázót. A maximális terhelést lásd a "Műszaki adatok" a 11. oldalon. A terhelés magában foglalja a rázógépre szerelt tálca, a tartozék alkatrészeknek és a mintáknak a súlyát.
- Mielőtt a rázót használatba veszi, győződjön meg arról, hogy megfelelően szerelték fel a tartozékokat. A "Tartozékok" a 88. oldalon utasításait be kell tartani.



VIGYÁZAT

Ujj- vagy kézsérülések veszélye leeső burkolat következtében.

 A rázógép burkolata véletlenül leeshet, és becsípheti vagy összepréselheti az ujjakat vagy a kezeket, ami sérülést okozhat. Ez akkor fordulhat elő, ha a burkolat rugója meghibásodik, és nem képes teljes mértékben nyitott helyzetben tartani a burkolatot. Ne működtesse a rázógépet hibás burkolatrugóval! Vegye fel a kapcsolatot a Thermo Fisher Scientific ügyfélszolgálatával, és haladéktalanul cseréltesse ki a burkolat rugóját.

Betöltés

Biztosítsa, hogy a töltet, a mintákat és tartozékokat is beleértve, a tálca közepéhez képest szimmetrikusan legyen elosztva. Kétlépcsős tálca használata a teljes kapacitás kihasználása nélkül: Optimális eredményt első sorban az alsó tálca feltöltésével és a közepéhez képest szimmetrikus elrendezésével lehet elérni.

Hőmérséklet-szabályozott rázógép használata esetén ügyeljen arra, hogy a rakományt úgy helyezze el, hogy a burkolat zárásakor és a burkolat zárt állapotában is elegendő hely maradjon az ütközés és a rakomány esetleges károsodásainak elkerüléséhez.



101. ábra: Példa a helyesen feltöltött tálcára

$\bigcirc \cdots \cdots \cdots \cdots \cdots \bigcirc \bigcirc$	n)
	12
·····································	Ŧ
	-
	• 🏚
· · · · · · · · · · · · · · · · · · ·	÷
	-
	•))))
$ \cdot \cdot$	Ŧ
	Ŧ
· · · · · · · · · · · · · · · · · · ·	•)))
[· · · · · · · · · · · · · · · · · · ·	Ŧ

102. ábra: Példa a hibásan feltöltött tálcára

MEGJEGYZÉS A maximális töltet tartalmazza a rázóra szerelt tálca, tartozékok és a minták súlyát. A maximális terhelésre vonatkozóan további információ az "Műszaki adatok" a 11. oldalon.

Rendeltetésszerű használat

A fordulatszám és a töltet határértékeire vonatkozó információk csak irányértékek a termék használatához. A csendes, kevés rezgéssel járó vagy rezgésmentes működés számos tényezőtől függ, például a töltési mennyiségtől, a lombiktartók és szőnyegek típusától és állapotától, az edény típusától, az edények elrendezésétől a tálcán, a rázógép alapjától és a névleges fordulatszámtól.

Mikor új vagy ismeretlen kombinációt állít be, vagy az alábbi paraméterek valamelyikét megváltoztatja: Lassan emelje 175 ford./perc fölé a fordulatszámot és ellenőrizze, hogy nem lép-e fel a töltet vagy a rázó nem kívánt lengése. Az üzemeltető felelős a rázó biztonságos üzemeltetéséért.

A Solaris rázógépeket kifejezetten alacsony hőleadásra és alacsony rezgésre tervezték, így ideálisak számos laboratóriumi berendezésben, például klímakamrákban, inkubátorokban és laboratóriumi hűtőszekrényekben történő használatra. A különböző speciális környezeti feltételek következtében, mint például a tálcák és a tartószerkezetek stabilitása, a felhasználó felelős a laboratóriumi berendezések biztonságos működéséért. Növelje meg lassan a fordulatszámot, és figyeljen a készülék szokatlan mozgására.

Biztonságos fordulatszám

A Solaris sorozatú rázógépek terheléstől független, biztonságos fordulatszáma az 56. táblázatban található. A töltet tartalmazza a rázóra szerelt tálca, tartozékok és a minták súlyát. A maximális terheléssel kapcsolatos további információkat lásd a "Műszaki adatok" a 11. oldalon.

Típus	Max. biztonságos fordulatszám
Solaris 2000	175 ford./perc
Solaris 4000	175 ford./perc
Solaris 2000 I	250 ford./perc
Solaris 2000 R	250 ford./perc
Solaris 4000 I	250 ford./perc
Solaris 4000 R	250 ford./perc

56. táblázat: Max. biztonságos fordulatszám

Maximális fordulatszám

A Solaris sorozat rázógépeit a legnagyobb, 525 fordulat/perc fordulatszámmal az 57. táblázat szerinti korlátozott terheléssel lehet csak üzemeltetni. A töltet tartalmazza a rázóra szerelt tálca, tartozékok és a minták súlyát. Vegye figyelembe az alkalmazott lombiktartók és edények megengedett fordulatszám tartományát, amik nem biztos, hogy kompatibilisek ezzel a fordulatszámmal.

Típus	Max. biztonságos terhelések
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

57. táblázat: Maximális terhelés maximális fordulatszámnál

MEGJEGYZÉS Mindig vegye figyelembe, hogy a növekvő fordulatszámnál a lombiktartók nyílni kezdenek és az edények fokozott mozgása révén a rendszer további átmeneti lengését okozzák.

Súly/fordulatszám jelleggörbék

Az alábbi súly/fordulatszám jelleggörbék segítséget nyújtanak a tervezett használathoz ideális fordulatszám/terhelés kombináció beállításához. A zöld tartomány azokat a fordulatszám - töltet kombinációkat mutatja, amik a rázás közben semmilyen vagy csak csekély lengést okoznak. A piros tartomány azokat a fordulatszám - töltet kombinációkat mutatja, amik a rázás közben semmilyen vagy csak csekély lengést okoznak. A piros tartomány azokat a fordulatszám - töltet kombinációkat mutatja, amik a rázás közben erős lengést és ezzel a rázó nem kívánt elmozdulását okozhatják. Mivel ezek a diagramok csak tájékoztatásul szolgálnak, vigyázzon amikor a fordulatszám - töltet kombináció közelíti a piros tartományt.

Vegye figyelembe, hogy a speciális alkalmazási körülményei a töltet és a készülék nem kívánt viselkedését okozhatják, mielőtt elérnének a berajzolt határvonalhoz. Lépésenként emelje a fordulatszámot, hogy láthassa speciális töltetének viselkedését. A töltet tartalmazza a rázóra szerelt tálca, tartozékok és a minták súlyát.

MEGJEGYZÉS Az alábbi ábrákon valamennyi kék vízszintes vonal egy adott tálca üres tömegét jelzi.



Solaris 2000

103. ábra: Solaris 2000 - Rendeltetésszerű használat

Solaris 4000



104. ábra: Solaris 4000 - Rendeltetésszerű használat



Solaris 2000 I / 2000 R

105. ábra: Solaris 2000 I / 2000 R - rendeltetésszerű használat

Feltöltés és rendeltetésszerű használat



106. ábra: Solaris 4000 I / 4000 R - rendeltetésszerű használat

4. Karbantartás és ápolás



FIGYELMEZ-TETÉS

Veszélyes anyagok kezelése okozta veszély

- Amennyiben veszélyes anyagot ráz, vegye figyelembe a "Laboratory Biosafety manual" kézikönyvet (a WHO kiadványa) és a rá vonatkozó nemzeti ajánlásokat. Ha a WHO "Laboratory Biosafety Manual" szerint II-es kockázati csoportba tartozó mikrobiológiai mintákat ráz, használjon aeroszoltömör biotömítéseket. A "Laboratory Biosafety Manual" a WHO internetes portálján érhető el (www.who.int). A magas rizikócsoportba tartozó anyagok esetén hozzon további biztonsági intézkedéseket.
- Ha a rázógépet vagy annak alkatrészeit toxinok vagy kórokozó anyagok szennyezték be, megfelelő fertőtlenítési intézkedéseket kell végrehajtania ("Dekontaminálás" a 124. oldalon; "Fertőtlenítés" a 124. oldalon).
- Veszélyhelyzet kialakulása esetén szakítsa meg a rázó áramellátását és azonnal hagyja el a területet.



FIGYELMEZ-

TETÉS

Fertőző anyagok okozta egészségkárosodás

Ha véletlenül folyadékok vagy más anyagok kerülnek a tálca alá, azonnal kapcsolja ki a rázógépet, válassza le a hálózatról és távolítsa el a tálcát (lásd "Tálcák" a 18. oldalon).

A szabványszerű laboratóriumi eljárásnak megfelelően mossa fel a kiömlött folyadékokat. Viseljen megfelelő személyi védőfelszerelést.

Az összes belső beállítást és javítást csak az arra feljogosított szerviztechnikussal végeztesse el. A felhasználó nem nyithatja ki a rázó házát.

Kövesse az adott tartozékokhoz mellékelt termékinformációkat, amik speciális tájékoztatást nyújtanak a szabályos tisztításról és karbantartásról. A jelen fejezet információit csak irányelvként használja.

4.1. Alapok

A személyek, anyagok és a környezet védelme érdekében köteles a rázót és tartozékait rendszeresen tisztítani és szükség esetén fertőtleníteni.

A Thermo Fisher Scientific javasolja, hogy legalább havonta egyszer tisztítsa meg és kézzel fertőtlenítse a laboratóriumi rázót. A helyiségek normál levegőjében ezerszám keringenek mikroorganizmusok, amik megtelepedhetnek a rázóban és veszélyeztethetik az Ön kultúráit.

A Thermo Fisher Scientific javasolja, hogy használjon 70%-os etanolt, 70%-os izo-propilalkoholt, 10%-os vagy hígabb ammónia oldat alapján.

MEGJEGYZÉS A behatoló folyadék károsíthatja a szerkezetet. Ne hagyja, hogy folyadékok, különösen is szerves oldószerek kerüljenek a szerkezetbe vagy annak csapágyához. A szerves oldószerek tönkre teszik a szerkezet csapágyának kenőzsírját. Ha folyadékok kerülnek a rázógép tisztítás céljából nem hozzáférhető részeibe, ne működtesse tovább a rázógépet, és lépjen kapcsolatba az ügyfélszolgálattal.

MEGJEGYZÉS A nem megengedett eljárások vagy anyagok megtámadhatják a rázó anyagait és hibás működéshez vezethetnek. Ne használjon más tisztítási vagy dekontaminációs eljárást, mint az itt leírt, ha nem biztos abban, hogy alkalmas az adott anyagokhoz. Csak olyan tisztítószereket használjon, amik nem károsítják az anyagokat. Kétség esetén forduljon a tisztítószer gyártójához.

MEGJEGYZÉS A hőmérséklet-szabályozott rázógépeknél nyitva kell hagyni a burkolatot, hogy el tudjon párologni a tisztítószer és a víz.

- Tisztítás, fertőtlenítés vagy dekontamináció előtt húzza ki a hálózati csatlakozót.
- Tisztítás, fertőtlenítés vagy dekontamináció előtt távolítsa el a rázóra szerelt tartozékokat és tálcá(ka)t.
- Az anyagok tisztításához használjon meleg vizet enyhe tisztítószerrel és puha ruhával. Kétség esetén forduljon a Thermo Fisher Scientifichez. Ez után tiszta vízzel öblítse le és alaposan szárítsa meg.
- Soha ne használjon maró tisztítószereket, pl. foszforsavat, hipót vagy súrolóport.
- Csak 6-8 közötti pH értékű fertőtlenítőszereket használjon.
- A kiömlött folyadékot haladéktalanul távolítsa el egy nem maró hatású tisztítószerrel megnedvesített szöszmentes ruhával (lásd a tisztítószer gyártójának utasításait).
- A kiömlött folyadék a tálca alá folyhat. Ha a tálca alatt kiömlik a folyadék, szerelje le a tálcát, és alaposan tisztítsa meg a tálca alatti tálat. Szükség esetén fertőtlenítse vagy mentesítse a szennyeződésektől.
- Ellenőrizze a rázó részegységeit és távolítsa el a kifolyt közeget és a szennyeződéseket.

A tartozékok ellenőrzése

MEGJEGYZÉS Ne használjon olyan rázót vagy tartozékot, amin sérülés jele látszik. Ajánlott rendszeres rutinellenőrzés során vizsgáltatni a tartozékokat, hogy garantálja a biztonságot.

A tartozékok alapos tisztítása után ellenőrizze azok sértetlenségét, elhasználtságát és korrózióját.

Fém részek

Ha károsodásra utaló jeleket, például korróziót, kopást vagy repedést észlel, haladéktalanul vonja ki a forgalomból a tartozék alkatrészeket.

Műanyag részek

Ellenőrizze a repedések jeleit, a kifakulásokat, horzsolásokat és töréseket. Meghibásodás jele esetén a vizsgált alkatrészt azonnal vonja ki a használatból.

Öntapadós szőnyegek

Minden használat előtt ellenőrizze az öntapadós szőnyegeket az elhasználódás vagy károsodás jelei, például repedések, elszíneződés, sárgulás, törékenység, alakváltozás, felületi kopás vagy kémiai hatások szempontjából. A kopás vagy károsodás jeleit mutató öntapadós szőnyegeket haladéktalanul ártalmatlanítsa.

4.2. Tisztítás

MEGJEGYZÉS A tisztítási folyamat megkezdése előtt a használó győződjön meg a tisztítószer gyártójánál arról, hogy tervezett folyamat nem károsítja az anyagokat.

MEGJEGYZÉS A behatoló folyadék károsíthatja a szerkezetet. Ne hagyja, hogy folyadékok, különösen is szerves oldószerek kerüljenek a szerkezetre vagy annak csapágyához. A szerves oldószerek tönkre teszik a szerkezet csapágyának kenőzsírját. Mossa le a készülék külsejét puha ruhával, illetve szappan és víz enyhe oldatával, öblítse le tiszta vízzel, majd alaposan szárítsa meg.

A rázógép és a használt tartozék alkatrészek megfelelő tisztítására vonatkozó további információkért lásd az "Alapok" a 122. oldalon.

MEGJEGYZÉS A hőmérséklet-szabályozott rázógépeknél nyitva kell hagyni a burkolatot, hogy el tudjon párologni a tisztítószer és a víz.

Kezelőfelület

A kezelőfelületet az alábbiak szerint tisztítsa meg:

- 1. Húzza ki a hálózati csatlakozót.
- 2. A kezelőfelületet száraz, mikroszálas kendővel törölje le.
- 3. Ha szükséges, nedvesítse meg a mikroszálas kendőt és ismét törölje át a kezelőfelületet.
- 4. Hagyjon minden részegységet megszáradni a levegőn.

4.3. Fertőtlenítés

FIGYELMEZTETÉS A veszélyes anyagok kezelése veszélyes. Ne érintse meg a fertőzött részeket. Súlyos fertőzést kaphat el a fertőzött részek megérintése esetén. A fertőző anyag az edény törése vagy kiöntés révén kerülhet a rázóba. Kontamináció esetén biztosítsa, hogy senki se legyen veszélyeztetve. Az érintett részeket azonnal fertőtlenítse.

MEGJEGYZÉS A nem megfelelő fertőtlenítő eljárás dekontamináló eljárás vagy szer károsíthatja az anyagokat. Győződjön meg arról, hogy a fertőtlenítési eljárás vagy szer nem árt az anyagoknak. Kétség esetén forduljon a fertőtlenítőszer gyártójához. Tartsa be az alkalmazott fertőtlenítőszer biztonsági és alkalmazási előírásait.

- Minden részegységet és felületet töröljön le 70%-os etanollal, a fertőtlenítés szükséges intenzitásától függően. Nem kerülhet nedvesség a szabadon lévő elektronikus alkatrészeket tartalmazó területekre!
- 2. Hagyjon minden részegységet megszáradni a levegőn.

MEGJEGYZÉS A hőmérséklet-szabályozott rázógépeknél nyitva kell hagyni a burkolatot, hogy el tudjon párologni a tisztítószer és a víz.

Saját maga felelős azért, hogy elérjék a követelményeinek megfelelő dekontaminációs fokot.

4.4. Dekontaminálás

FIGYELMEZTETÉS A veszélyes anyagok kezelése veszélyes. Ne érintse meg a kontaminált részeket. A szennyezett alkatrészek megérintésekor szennyeződés történhet. A kontaminált anyag az edény törése vagy kiöntés révén kerülhet a rázóba. Kontamináció esetén biztosítsa, hogy senki se legyen veszélyeztetve. Haladéktalanul dekontaminálja az érintett részeket.

MEGJEGYZÉS A nem megfelelő dekontamináló eljárás vagy szer károsíthatja az anyagokat. Győződjön meg arról, hogy a szennyeződésmentesítési módszer vagy szer nem károsítja az anyagokat. Kétség esetén forduljon a dekontamináló szer gyártójához. Tartsa be az alkalmazott dekontamináló szer biztonsági és alkalmazási előírásait.

A Thermo Fischer Scientific az alábbi módszert ajánlja.

- Minden részegységet és felületet töröljön le 70%-os etanollal. Ne nedvesítse meg a szabadon levő elektronikus részeket.
- 2. Hagyjon minden részegységet megszáradni a levegőn.

MEGJEGYZÉS A hőmérséklet-szabályozott rázógépeknél nyitva kell hagyni a burkolatot, hogy el tudjon párologni a tisztítószer és a víz.

Saját maga felelős azért, hogy elérjék a követelményeinek megfelelő dekontaminációs fokot.

4.5. Autoklávozás

MEGJEGYZÉS Soha ne lépje túl az autoklávozási hőmérséklet és időtartam megengedett értékeit! Nem megengedett a vegyi adalékok használata a gőzben.

Az autoklávozás előtt:

- Szereljen le minden tartozékot.
- Alaposan öblítse le a vegyszerek vagy fertőtlenítőszerek minden maradványát. Végül desztillált vízzel öblítse át.

További információkat az autoklávozhatóságról az alábbi táblázatban talál:

Cikk	Autoklávozható	Műszaki adatok
Rázó	Nem	-
Tálca	Igen	121 °C, 20 perc
Lombiktartó (rugós hüvelyek nélkül)	Igen	121 °C, 20 perc
Csavarhúzó	Nem	-
Lombiktartó Mikrotiter/Deepwell tálcákhoz	lgen	121 °C, 20 perc
Tapadószőnyeg	Nem	-
Főzőpohár állvány	lgen	121 °C, 20 perc
Szétválasztó tölcsértartó	lgen	121 °C, 20 perc
Kémcsőállvány – fém alkatrészek	Igen	121 °C, 20 perc
Kémcsőállvány – műanyag alkatrészek	Nem	-
Kémcsőállvány – habbetétek	Nem	-
Állítható edénytartó – fém alkatrészek	lgen	121 °C, 20 perc
Állítható edénytartó – csúszásgátló alátét	Nem	-
Többfunkciós tároló – fémtároló	lgen	121 °C, 20 perc
Többfunkciós tároló – gumiszőnyeg	Nem	-
Állítható szögű kémcsőállvány tartó – fém alkatrészek	lgen	121 °C, 20 perc
Állítható szögű kémcsőállvány tartó – műanyag alkatrészek	Nem	-
Állítható szögű kémcsőállvány tartó – habbetétek	Nem	-

58. táblázat: Az anyagok autoklávozhatósága

Győződjön meg arról, hogy a saját követelményeinek megfelelő sterilitást elérték-e.

Az autoklávozás következtében kissé megváltozhat a megjelenés és a szín.

4.6. Hőmérsékletkalibrálás

A hőmérséklet-szabályozott rázógép hőmérsékletszabályozási pontja egy hőmérsékletmérővel a szükséges mérési pontosságig kalibrálható, hogy garantálja az inkubált (csak fűtött) vagy hűtött (hűtött és fűtött) rázási folyamat teljes reprodukálhatóságát.

4. 6. 1. Bevált hőmérsékletkalibrálási eljárások

A rázógép hőmérsékletkalibrálása előtt és alatt gondosan tartsa be az utasításokat. Bármilyen zavar, vagy az ajánlott legjobb gyakorlattól való eltérés azt okozhatja, hogy a minta hőmérséklete nem stabilizálódik, és a kalibrálás sikertelen lesz.

A bevált gyakorlatok a következők:

- Használjon 250 ml űrtartalmú lombikot.
- Töltse meg a lombikot pontosan 100 ml vízzel.
- Zárja le teljesen a burkolatot, és tartsa zárva a kalibrálás egésze alatt.
- Hagyjon elegendő időt annak garantálásához, hogy a minta hőmérséklete teljes mértékben stabilizálódjon.
- Győződjön meg arról, hogy a környezeti feltételek, mint például a hőmérséklet és a páratartalom változatlanok maradnak a kalibrálás alatt. Már apró változások, például egy ajtó vagy ablak kinyitása vagy a légkondicionáló beállításainak megváltoztatása is befolyásolhatja a folyamatban lévő kalibrációs folyamat eredményét.
- A rázógépet évente egyszer ajánlott kalibrálni.

4. 6. 2. A hőmérsékletkalibrálás végrehajtása

A kalibráláshoz a következőképpen járjon el:

- 1. A kalibrálás háromféleképpen indítható el:
 - » Az "Alap beállítás" a 54. oldalon leírt alapbeállítási eljárás végén lehetősége van arra, hogy az alapbeállítási eljárásból történő kilépés helyett elvégezze a hőmérsékletkalibrálást.
 - » A rázógép megjeleníti az éves kalibrálásra vonatkozó emlékeztetőt.
 - » Ezután a navigációs sáv Beállítások ikonjának, majd a megjelenő Hőmérsékletkalibrálás képernyő alsó részén található Kalibrálás gombnak a megérintésével indíthatja el a kalibrálást.
 - » A navigációs sáv Beállítások ikonjának, majd a Kalibrálás gombnak a megérintésével szükség esetén elindíthatja a folyamatot.

MEGJEGYZÉS A kalibrálási emlékeztető kikapcsolható a Kalibrálási üzenetek kikapcsolása jelölőnégyzet bejelölésével a Beállítások oldalon. 2. Várjon, amíg megjelenik a Hőmérsékletkalibrálás képernyőn látható 107. ábra.



107. ábra: Kalibrálás: Kezdőképernyő

3. A folyamat elindításához érintse meg a Kalibrálás elindítása gombot.



108. ábra: Kalibrálás: Erlenmeyer-lombik feltöltésének kérése

- Kövesse a képernyőn megjelenő utasításokat: Töltsön meg egy 250 ml-es Erlenmeyer-lombikot 100 ml vízzel.
- Helyezze a lombikot egy lombiktartóba, és rögzítse a lombiktartót körülbelül a rázótálca közepén.
- Akassza a lombikba a kívánt mérési pontossággal rendelkező hőmérsékletmérő eszköz szondáját. A szondának a folyadékba kell merülnie, de nem szabad hozzáérnie a lombik aljához vagy oldalához.
- 7. Zárja le a lombikot, hogy elkerülje a folyadék kiömlését.
- 8. Zárja le a burkolatot a rázótálca fölött.

9. A folytatáshoz érintse meg a Tovább gombot.



109. ábra: Kalibrálás: Kalibrálás megkezdésére irányuló felszólítás

MEGJEGYZÉS A rázógép a Start gomb megérintésekor azonnal elkezdi a rázást.

10. Kövesse a képernyőn megjelenő utasításokat: Érintse meg a Start gombot az előre beállított hőmérséklet használatához, vagy érintse meg a Hőmérséklet névleges értéke mezőt, és adjon meg egy másik kívánt hőmérsékletet, mielőtt megérinti a Start gombot.

A képernyőn látható, hogy miként történik a rázókamra melegítése (vagy hűtése) a kiválasztott kalibrációs hőmérsékletre.



110. ábra: Kalibrálás: A kamra kalibrációs hőmérséklet szerinti stabilizálása

 Várjon legalább két és fél órát (150 perc), hogy a kamra elérje a hőmérsékletmérés szempontjából stabil állapotot.

A stabil állapotot az **A mintakamra hőmérséklete stabilizálódott** üzenet jelzi, amint azt az alábbi 111. ábra mutatja.



111. ábra: Kalibrálás: A kamra kalibrációs hőmérséklet szerinti stabilizálása sikeresen megtörtént

- 12. Kövesse a képernyőn megjelenő utasításokat:
 - » Amint megjelenik a 111. ábran látható képernyő, rendszeresen hasonlítsa össze a hőelem hőmérsékletértékét. Ha a hőmérsékletet stabilnak találja, a Szerkesztés gomb megérintésével korrigálhatja, vagy a Befejezés gombot megérintve befejezheti a kalibrálási folyamatot.



112. ábra: Kalibrálás: A rázógép hőmérséklet-kijelzésének korrigálása

- A megjelenő beállítókerék segítségével korrigálja a rázógép által kijelzett hőmérsékletet a mérőműszer szerinti értékre.
- 14. A módosítások mentéséhez érintse meg a Mentés gombot.

15. A Hőmérsékletkalibrálás befejeződött üzenet azt jelzi, hogy a folyamat sikeresen befejeződött.

<	Shake Lab 1 11.22.24AM Temp Calibration	
۵.	Step 5 of 5	
首	Temperature calibration complete.	
Ē		
	Re-Calibrate	()

113. ábra: Kalibrálás: A hőmérsékletkalibrálás sikeresen befejeződött

A kalibrálási folyamat megismétléséhez érintse meg az Újrakalibrálás gombot.

4. 6. 3. A hőmérsékletkalibrálás lekérdezése

Az aktuális kalibrációs hőmérséklet bármikor előhívható:

- Érintse meg a navigációs sáv Beállítások ikonját, majd a megjelenő Hőmérsékletkalibrálás képernyő alsó részén található Kalibrálás gombot.
- Megjelenik a 107. ábran látható Kalibrálás képernyő. Ez a képernyő az aktuális Kalibrált rázóhőmérsékletet mutatja.

MEGJEGYZÉS Hőmérséklet-szabályozott rázógépek esetében a kalibrációs hőmérsékletet arról a képernyőről is leolvashatja, ahol a hőmérsékletet beállította. Egy példa található a "A hőmérséklet beállítása" a 64. oldalon.

4. 7. A készülékszoftver telepítése

Ha a Fájlok és információk képernyőn megérinti a Készülékszoftver gombot (lásd a "Fájlok és info" a 77. oldalon), a rendszer felszólítja a rendszergazda szintű hozzáférési kód megadására, és megjelenik a 114. ábra szerinti készülékszoftver képernyő. Ezen a képernyőn keresztül új készülékszoftvert telepíthet USB-meghajtóról.

A szervizszerződésétől függően előfordulhat, hogy a helyszíni szerviztechnikustól kap egy letöltési hivatkozást az új készülékszoftverhez. Ügyeljen arra, hogy csak engedélyezett forrásból származó készülékszoftvert telepítsen. A nem engedélyezett forrásból származó szoftverek érvényteleníthetik a Thermofisher Scientific által vállalt garanciát.

- 1. Töltse le a ZIP fájlt a helyszíni szerviztechnikusától kapott hivatkozáson keresztül.
- 2. Csomagolja ki a ZIP fájlt egy üres USB-meghajtóra.

MEGJEGYZÉS Győződjön meg arról, hogy elegendő szabad hely áll rendelkezésre az USB-meghajtón a ZIP fájl kicsomagolásához.

3. Ha a rázógép továbbra is működik, állítsa le a kijelzőn látható Stop gomb megérintésével.

 Nyugtázza az esetlegesen fennálló riasztási üzeneteket az "Üzemállapot" a 66. oldalon leírtak szerint.

MEGJEGYZÉS Ha a frissítést a rázógép működése közben kísérli meg elvégezni, megjelenik a következő üzenet: A frissítést nem szabad elvégezni, amíg a rázógép rázó üzemmód szerint működik. Ez az üzenet tartalmaz egy Rázó leállítása gombot, amely lehetővé teszi a futás leállítását és a frissítést.

- 5. Nyomja meg a navigációs sávon a Fájlok és infó jelet.
- A Fájlok és információk képernyőn görgessen egészen le a Készülékszoftver gomb megjelenítéséhez.
- Érintse meg a Készülékszoftver gombot a Készülékszoftver képernyő megnyitásához, amely az aktuálisan telepített készülékszoftver-verziókat jeleníti meg.
- Csatlakoztassa hozzá az USB-meghajtót a rázógép USB-csatlakozóinak egyikéhez. Az USBcsatlakozók helyei a "Termék áttekintés" a 44. oldalon.
- 9. Adja meg a rendszergazda szintű hozzáférési kódot.



114. ábra: Készülékszoftver képernyő

10. Érintse meg a Készülékszoftver-frissítés végrehajtása gombot.

MEGJEGYZÉS Ha a rázógép nem talál USB-meghajtót, megjelenik a következő kérés: Kérjük, a frissítés elindításához helyezzen be egy USB-meghajtót a telepítendő készülékszoftverrel. Ekkor helyezze be az USB-meghajtót, vagy ellenőrizze, hogy megfelelően csatlakoztatta-e hozzá az USB-csatlakozóhoz, és javítsa ki a problémát.

11. Készülékszoftver-frissítés elérhetővé válása esetén megjelenik a készülékszoftver-frissítés képernyőn a 115. ábra szerinti lista. Érintse meg a frissíteni kívánt készülékszoftver elemet.

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

115. ábra: A telepítendő készülékszoftver elem kiválasztása

MEGJEGYZÉS A legfrissebb verziójú készülékszoftver elemek szürkén jelennek meg, és nem választhatók ki. Ha nincsenek frissítendő elemek, akkor a **Jelenleg nincs új frissítés a szoftverhez** üzenet jelenik meg.

- 12. Érintse meg a Kiválasztás gombot.
- Ha a memóriakártya egynél több készülékszoftver-verziót tartalmaz, válassza ki a megfelelőt. Érintse meg a kívánt készülékszoftver-verziót, majd pedig a Kiválasztás gombot.



116. ábra: A telepítendő készülékszoftver-verzió kiválasztása

MEGJEGYZÉS Ha nem biztos benne, hogy melyik verziót kell telepítenie, kérjük, forduljon a letöltési hivatkozást biztosító technikushoz. Megjelenik azoknak az elemeknek a listája, amelyeknél frissítést kell végrehajtani. A frissítés elindításához érintse meg az OK gombot.

Firmware Upgrade	>
Current Version: 176	Upgrade Version: 254
require a rebo	ot of the interface.

117. ábra: A frissítendő elemek listája

MEGJEGYZÉS A felhasználói felület kezeli Ön helyett a kölcsönös függőségeket: Kiválasztja a frissítendő készülékszoftver elemet, és a szoftver automatikusan hozzáadja a telepítéshez szükséges összes elemet.

MEGJEGYZÉS Ha a 115. ábra listájából a H.M.I. lehetőséget csak a grafikus felhasználói felület készülékszoftverének frissítése céljából jelöli ki, előfordulhat, hogy a grafikus felhasználói felület kijelzője a készülékszoftver-frissítés megkezdése után akár egy percen keresztül sem reagál a felhasználói felület aktív területének megérintésére. Kérjük, hogy ez idő alatt <u>ne</u> kapcsolja ki a rázógépet, ellenkező esetben a frissítés nem fejeződik be sikeresen. A rázógép egy rövid idő elteltével újraindul, majd ismét kész a használatra.

MEGJEGYZÉS Ha két vagy három lehetőséget választ ki a 115. ábran látható listából <u>több</u> készülékszoftverösszetevő frissítéséhez, a rázógépet a frissítési folyamat közben semmiképpen sem szabad kikapcsolni.

 Különböző üzenetek jelennek meg, amelyek a telepítés előrehaladásáról tájékoztatják Önt a 118. ábran látható módon.



118. ábra: Készülékszoftver-frissítés telepítésével kapcsolatos üzenetek

 Ha megjelenik a rázógép újraindítására vonatkozó felszólítás, kapcsolja ki, majd újra be a készüléket a hálózati kapcsolóval.

4.8. Tálcacsere

A tálca összeszereléséhez szükséges alapvető tudnivalókat lásd a "A tálca felszerelése" a 89. oldalon. Az egyes rázógépekhez tartozó cseretálcák teljes listáját lásd az "Tálcák" a 18. oldalon.

MEGJEGYZÉS A hőmérséklet-szabályozott rázógépeket kizárólag a gyárilag előre beszerelt tálcákkal szabad üzemeltetni.

4.9. Szerviz

A Thermo Fischer Scientific ajánlja, hogy évente egyszer végeztessen karbantartást a rázón és tartozékain egy feljogosított szerviztechnikussal. A szerviztechnikus az alábbiakat ellenőrzi:

- Elektromos felszerelések
- A telepítési hely alkalmassága
- Biztonsági rendszer
- Alkalmazott tartozékok
- A lombiktartó, tálcák és a rázó egyéb tartozékai rögzítése

A szerviz előtt alaposan meg kell tisztítani és dekontaminálni a rázót, hogy biztosítsa a teljes és biztonságos belső felülvizsgálatot.

Erre a szoilgáltatásra a Thermo Fischer Scientific felülvizsgálati és karbantartási szerződést kínál. Az esetleg szükséges javításokat a garancia keretében díjtalanul, a garanciális időn túl díj ellenében végezzük el. Ez csak akkor érvényes, ha a készülékbe csak a Thermo Fischer Scientific szerviztechnikusai avatkoztak be.

4.10. Elküldés és megsemmisítés

FIGYELMEZTETÉS Fertőző anyagok okozta egészségkárosodás. Ha a rázót és tartozékait használaton kívül kell helyezze és megsemmisítse, meg kell tisztítsa és ha szükséges fertőtlenítse vagy dekontaminálja. Kétség esetén forduljon a Thermo Fischer Scientific ügyfélszolgálatához.

A rázó megsemmisítése során tartsa be országának előírásait. A rázó megsemmisítése ügyében forduljon a Thermo Fischer Scientific ügyfélszolgálatához. Az elérhetőségi információkat ennek az útmutatónak a hátoldalán találja, vagy az interneten érheti el a <u>http://thermofisher.com</u> címen.

Lásd a szállításra és szállítmányozásra vonatkozó információkat ("Szállítás" a 52. oldalon és "Elküldés" a 59. oldalon).

5. A hibák okai és elhárításuk

Sz.	Megnevezés	Megoldások	Jel
Hiba			
11999	Belső hiba lépett fel.	Belső hiba lépett fel. Kérjük, indítsa újra a készüléket a hálózati kapcsoló ki- és bekapcsolásával. Ha a hibajelzés továbbra is megjelenik, forduljon egy szerviztechnikushoz.	
Figyelmezt	etések		
3	Hőmérsékletkalibrálás esedékes. Eltelt 1 év.	Az egyéves érvényességi idő lejárt. Ismételje meg a hőmérsékletkalibrálást a "Hőmérsékletkalibrálás" a 126. oldalon leírtak szerint.	CAL
4	Áramellátás helyreállítva, az automatikus újraindítás bekapcsolva.	A legutóbbi futás során áramkimaradás volt. Az áramellátás helyreállítása után a működés automatikusan folytatódott.	Ę
6	A környezeti hőmérséklet érzékelői meghibásodtak.	A környezeti hőmérséklet érzékelői által mért hőmérséklet valószínűtlen.	
10	Az 1. ventilátor fordulatszáma meghaladja a határértéket.	Az 1. ventilátor fordulatszáma a valószínűségi határérték felett van.	•2
11	Az 2. ventilátor fordulatszáma meghaladja a határértéket.	Az 2. ventilátor fordulatszáma a valószínűségi határérték felett van.	•2
12	Az 3. ventilátor fordulatszáma meghaladja a határértéket.	Az 3. ventilátor fordulatszáma a valószínűségi határérték felett van.	•2
13	Az 4. ventilátor fordulatszáma meghaladja a határértéket.	Az 4. ventilátor fordulatszáma a valószínűségi határérték felett van.	•2
20	A burkolat már régóta nyitva van.	A kamra hőmérsékletének fenntartása érdekében zárja le a rázógép burkolatát.	4
Riasztások			
1	Túlmelegedési riasztás.	A kamra hőmérséklete túllépte a felső riasztási értéket. Kérjük, ellenőrizze a mintákat, a környezeti feltételeket, és/vagy módosítsa a beállításokat.	
2	Túl alacsony hőmérsékletre figyelmeztető riasztás.	A kamra hőmérséklete túllépte az alsó riasztási értéket. Kérjük, ellenőrizze a mintákat, a környezeti feltételeket, és/vagy módosítsa a beállításokat.	
3	Sikertelen volt az áramkimaradás utáni automatikus újraindulás.	A legutóbbi futás során áramkimaradás volt. Nem sikerült az automatikus újraindítás végrehajtása.	Ċ
5	Túl lassú a meghajtás gyorsulása. Nem érhető el a kívánt névleges fordulatszám.	Nem volt idejében elérhető a kívánt névleges fordulatszám. Ellenőrizze a beállításait és/vagy a tálcán levő töltetét (tartozékok és minták).	\odot
7	Az 1. ventilátor fordulatszámmérése váratlan leállást jelez.	Kapcsolja ki, majd ismét be a készüléket. Ha a hibajelzés továbbra is megjelenik, forduljon egy szerviztechnikushoz.	•2

Sz.	Megnevezés	Megoldások	Jel
8	Az 2. ventilátor fordulatszámmérése váratlan leállást jelez.	Kapcsolja ki, majd ismét be a készüléket. Ha a hibajelzés továbbra is megjelenik, forduljon egy szerviztechnikushoz.	3
9	Az 3. ventilátor fordulatszámmérése váratlan leállást jelez.	Kapcsolja ki, majd ismét be a készüléket. Ha a hibajelzés továbbra is megjelenik, forduljon egy szerviztechnikushoz.	3
10	Az 4. ventilátor fordulatszámmérése váratlan leállást jelez.	Kapcsolja ki, majd ismét be a készüléket. Ha a hibajelzés továbbra is megjelenik, forduljon egy szervíztechnikushoz.	3
12	Hiba a meghajtás indításánál - nem érkezett fordulatszámjel.	A rázó tálcája blokkolódott. Ellenőrizze, hogy a tálca körül elegendő hely van-e és/vagy csökkentse a tálcán levő töltetet (tartozékok és minták). Végül nyomja meg ismét a rázó START gombját. Ha a hibajelzés továbbra is megjelenik, forduljon egy szerviztechnikushoz.	8
24	Fordulatszámmérési hiba a működés közben.	Szokatlan fordulatszám-változást érzékelt. Ellenőrizze a töltetet (minták és tartozékok) és/vagy a lombiktartót a rázó tálcáján. Végül nyomja meg ismét a rázó START gombját. Ha a hibajelzés továbbra is megjelenik, forduljon egy szerviztechnikushoz.	4
26	A fordulatszámmérés működés közben nem várt leállást érzékelt.	Ellenőrizze, hogy a tálca körül elegendő hely van-e és/vagy csökkentse a tálcán levő töltetet (tartozékok és minták). Győződjön meg arról, hogy a rázogép hátulján található biztosíték nem oldott-e ki ("Biztosíték" a 47. oldalon). Végül nyomja meg ismét a rázó START gombját. Ha a hibajelzés továbbra is megjelenik, forduljon egy szerviztechnikushoz.	٢
82	A motor áramnak mérése túlterhelést érzékelt.	Motor túláramot érzékelt. Ne töltse fel vagy ürítse ki a tálcát a készülék működése közben. Csökkentse a fordulatszámot vagy illessze a tálcán levő töltetet (tartozékok és minták).	0
83	A motor áramának mért értékei a megengedett tartományon kívül vannak.	Motor túláramot érzékelt. Ne töltse fel vagy ürítse ki a tálcát a készülék működése közben. Csökkentse a fordulatszámot vagy illessze a tálcán levő töltetet (tartozékok és minták).	

59. táblázat: Hiba-, figyelmeztető és riasztási üzenetek listája

MEGJEGYZÉS Ha olyan hibajelzés jelenik meg, ami nem szerepel ebben a táblázatban, forduljon egy szerviztechnikushoz.

GPL (General Public License, nyilvános licenc)

A készülék szoftverének egy része GPL, LGPL vagy más nyílt forráskódú licenc alatt közzétett nyílt forráskódú szoftvereket használ. A táblázatban részletesen felsorolt könyvtárakról van szó. A felhasznált könyvtárak (harmadik féltől származó könyvtárak) forráskódja beszerezhető a Thermo Fisher Scientific cégtől, ha a megfelelő licenc ezt előírja. A felhasznált nyílt forráskódú szoftverek licencfeltételei a rendelkezésre bocsátott forráskódcsomag részét képezik.

Könyvtár	Verzió	Felhasználó	Licenc
Qt	5.8	BSP	LGPLv3
Log4Cplus	1.2.0	Alkalmazás	Apache Public Licence v2/két klauzulás BSD licenc
boost	1.72.0	Alkalmazás	Boost licenc 1.0
json11	1.0.0	Alkalmazás	MIT licenc

Harmadik féltől származó könyvtárak

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

Tárgymutató

Α

A dugattyútartó és az edény behelyezése 92 A főzőpohár állványok felszerelése 101 A hibák okai és elhárításuk 135 A kémcsőállvány felszerelése 96 A kezelési útmutatóban használt jelek 7 Alap beállítás 54 Állítható edénytartó felszerelése 99 Állítható szögű kémcsőállvány-tartó felszerelése 97 Alvó mód 75 Ápolás 121 A szállítmány tartalma 49 A szétválasztó tölcsértartó felszerelése 102 A tálca felszerelése 89 A tartozékok ellenőrzése 123 Autoklávozás 125 Automatikus újraindítás 73 Az öntapadós szőnyeg lerakása és használata 106

В

Beállítások 68 Betöltés 116 Biztonságos fordulatszám 117 Biztosíték 47

С

Csatlakozások 46

D

Dekontaminálás 124

Ε

Elküldés 59, 134 Eseménynapló 84 Ethernet 47

F

Fájlok és info 77 Feltöltés 115 Fém részek 123 Fényerő 73 Fertőtlenítés 124 Figyelmeztetések 69

G

Gázelosztó felszerelése 113 GPL (General Public License, nyilvános licenc) 137 Grafikonok 85 Grafikus felhasználói felület 60

Η

Hálózati csatlakozás 47 Hangerő 69 Hőmérsékletkalibrálás 126 Hőmérséklet névleges értéke 73

ldő 75 Időmód 73 Irányelvek 43

J

Jelek és kifejezések 6

Κ

Karbantartás 121 Készülék megnevezés 76 Kétlépcsős tálcák 91 Kezelés 60 Kijelző 73

L

Lombiktartó felszerelése Mikrotiter/Deepwell tálcákhoz 98

М

Maximális fordulatszám 117 Megsemmisítés 134 Műanyag részek 123 Műszaki adatok 11

Ν

Némítás időtartama 70 Névleges fordulatszám 73 Nyelv 74

Ρ

Programok 78

R

Régió 76 Rendeltetésszerű használat 5, 115, 117 Riasztások 69

S

Szabványok 43 Szállítás 52 Szögletes lombiktartó felszerelése 94

Т

Tálcák 18 Tartozékok 17, 88 Telepítési hely 50 Termék áttekintés 44 Tisztítás 123 Többfunkciós tároló felszerelése 105

U

Univerzális tálcák 90 USB 47 Üzemi paraméterek 73

thermo scientific



Gyártva a



Thermo Electron LED GmbH Zweigniederlassung Osterode Am Kalkberg, 37520 Osterode am Harz Germany Származási ország: 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 az eredeti kezelési útmutató. Ez a kezelési útmutató az eredeti Kezelési útmutató fordítása.

thermofisher.com/shaker

© 2019–2023 Thermo Fisher Scientific Inc. Minden jog fenntartva.

Ha nem jelöljük kifejezetten másképp, akkor a márkanevek a Thermo Fisher Scientific Inc. és a hozzá kapcsolódó cégek tulajdonát képezik. Nem minden termék kapható minden országban. Pontosabb információkért érdeklődjön helyi kereskedelmi partnerünknél.

A jelen kezelési útmutatóban közölt képet tájékoztató jellegűek. Az ott bemutatott beállítások és nyelvek eltérhetnek.

Ausztrália +61 39757 4300

Ausztria +43 1 801 40 0

Belgium +32 53 73 42 41

Kína +800 810 5118 vagy +400 650 5118

Franciaország +33 2 2803 2180

Németországban belföldről, díjtalan 0800 1 536 376

Németország külföldről +49 6184 90 6000 India +91 22 6716 2200

Olaszország +39 02 95059 552

Japán +81 3 5826 1616

Hollandia +31 76 579 55 55

Újzéland +64 9 980 6700

Északi országok/Baltikum/ FÁK országok +358 10 329 2200

Oroszország +7 812 703 42 15 Spanyolország/Portugália +34 93 223 09 18

Svájc +41 44 454 12 12

Nagy-Britannia / Írország +44 870 609 9203

USA/Kanada +1 866 984 3766

Egyéb ázsiai országok +852 2885 4613

Más országok +49 6184 90 6000

