

Thermo Scientific Míchadla konstrukční řady Solaris

Návody

70900203-d • 2023-03



Obsah

Úvod	5
O tomto návodu	
Použití v souladu s určením	
Otevřené míchadlo	
Inkubované míchadlo (jen vy	hřívané) a chlazené míchadlo
(chlazené a vyhřívané)	
Potřebné znalosti	
Signální slova a symboly	
Symboly na míchadle a příslu	ušenství
Symboly použité v návodu .	
Bezpečnostní pokyny	
1. Technické specifikace	
	44
Michaelo Thermo Scientific S	Joiaris 2000 12 Solaris 4000 12
Michadio Thermo Scientific S	
Michadio Thermo Scientific S	Jolaris 2000 I 13 Solaris 4000 I 14
Michaele Thermo Scientific S	Solaris 2000 D
Michaelo Thermo Scientific S	Solaris 2000 R
1. 2. 1. Zakladny	
1. 2. 2. Drzaky banek	
1. 2. 3. Drzaky banek pro mikrotitrachi.	/Deepweil desticky 23
1. 2. 4. Stojany na zkumavky	
1. 2. 5. Drzaky stojanu na zkumavky s	
1. 2. 6. Nastavitelne drzaky nadob	
1. 2. 7. Civercove drzaky hadob	
1. 2. 0. Stojany pro kadinky inalgene .	
1. 2. 9. DIZAKY PIO UCIICI HALEVKY	۵۵ ۵۵
1. 2. 11. Aufiezi il poulozky	40 ،

1. 3.	Normy a směrnice	43
1. 4.	Přehled výrobku	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. Přípojky	46
2. Přer	prava a instalace	49
2.1	Vybalení	49
	Obsah dodávky	49
2 2	Místo instalaço	50
2.2. 0.0		50
Z. J.	Maninulago e míchadlam	52
.		52
Z. 4.		53
2. 5.	Siťový konektor	53
2. 6.	Základní nastavení	54
2.7.	Skladování	59
2. 8.	Odeslání	59
3. Obs	luha	60
3 1	Zannutí/wynnutí	60
J. I. 2 0		00
3. Z.		0U
ა. ვ	2. 1. Nastavení nejoulezitejsích provozních parametru michadia	66
ט. ז	2. 2. FIOVOZIII Stav	68
3.	2.4 Programy	78
2 2	Dříclušanetví	88
3	3 1 Montáž základny	89
3	3 2 Nasazení držáků baněk a nádob	92
3.	3. 3. Instalace čtvercového držáku baněk	94
3.	3. 4. Instalace stojanu na zkumavky	96
3.	3. 5. Instalace držáku stojanu na zkumavky s nastavitelným úhlem	97
3.	3. 6. Instalace držáku pro mikrotitrační/DeepWell destičky	98
3.	3. 7. Instalace nastavitelného držáku nádob	99

3. 3. 8. Instalace stojanu pro kádinky	101
3. 3. 9. Instalace držáku pro dělicí nálevku	102
3. 3. 10. Instalace univerzálního odkládacího tácu	105
3. 3. 11. Položení a použití adhezní podložky	106
3. 3. 12. Instalace rozdělovače plynu	. 113
3. 4. Nakládání a použití v souladu s určením	115
Vkládání	. 116
Použití v souladu s určením	. 117
4. Údržba a péče	121
4. 1. Základní principy	122
Kontrola příslušenství	123
4. 2. Čištění	123
Ovládací panel	123
4. 3. Dezinfekce	124
4. 4. Dekontaminace	124
4. 5. Autoklávování	125
4. 6. Kalibrace teploty	126
4. 6. 1. Osvědčené postupy kalibrace teploty	126
4. 6. 2. Provedení kalibrace teploty	126
4. 6. 3. Vyhledání kalibrační teploty	130
4. 7. Instalace firmwaru	130
4. 8. Výměna základny	134
4. 9. Servis	134
4. 10. Odeslání a likvidace	134
5. Příčiny a odstraňování poruch	135
GPL (General Public License, obecná veřejná licence)	137
Rejstřík	138

Úvod

O tomto návodu

Tento návod k použití popisuje následující míchadla a jejich příslušenství ("1. 2. Příslušenství" na straně 17):

Míchadlo	Č. výrobku
Michadlo Thermo Scientific Solaris 2000, 100–240 V ± 10 %, 50 / 60 Hz	SK2000
Michadlo Thermo Scientific Solaris 4000, 100–240 V \pm 10 %, 50 / 60 Hz	SK4000
Michadlo Thermo Scientific Solaris 2000 I, 100–120, 200–240 V ±10 %, 50 / 60 Hz	SK2001
Michadlo Thermo Scientific Solaris 4000 I, 100–120, 200–240 V ±10 %, 50 / 60 Hz	SK4001
Michadlo Thermo Scientific Solaris 2000 R, 100–240 V ±10 %, 50 / 60 Hz	SK2002
Michadlo Thermo Scientific Solaris 4000 R, 100–240 V ±10 %, 50 / 60 Hz	SK4002

Popisy a pokyny v tomto návodu se nevztahují na úplný název výrobku, ale na jednotlivá míchadla podle tříd:

- Otevřené míchadlo
- Inkubované míchadlo (jen vyhřívané)
- Chlazené míchadlo (chlazené a vyhřívané)

Použití v souladu s určením

Otevřené míchadlo

Toto zařízení se používá ve výzkumu a výrobě pro míchání roztoků po dvourozměrné kruhové dráze, zároveň umožňuje regulovat aplikační otáčky a délku míchání.

Inkubované míchadlo (jen vyhřívané) a chlazené míchadlo (chlazené

a vyhřívané)

Toto zařízení se používá ve výzkumu a výrobě pro míchání roztoků po dvourozměrné kruhové dráze, zároveň umožňuje regulovat aplikační otáčky, délku míchání a teplotu.

Potřebné znalosti

Míchadlo smí provozovat jen proškolený personál.

Signální slova a symboly

Signální slovo	Stupeň nebezpečí
VÝSTRAHA	Upozorňuje na nebezpečnou situaci, která může vést k těžkým nebo smrtelným zraněním osob, jestliže nezabráníte jejímu vzniku.
POZOR	Upozorňuje na nebezpečnou situaci, která může vést k drobným nebo lehkým zraněním osob, jestliže nezabráníte jejímu vzniku.
UPOZORNĚNÍ	Upozorňuje na důležité informace, které se nevztahují k nebezpečím.

Symboly na míchadle a příslušenství

Dodržte pokyny v návodu, aby nedošlo k ohrožení vás ani okolí.

8	Další informace najdete v návodu k použití
	Vytáhněte zástrčku z elektrické sítě
	Obecné nebezpečí
	Nebezpečí popálení na horkém povrchu
	Nebezpečí zranění prstů nebo rukou při neúmyslném zavření mechanických dílů

Symboly použité v návodu

Dodržte pokyny v návodu, aby nedošlo k ohrožení vás ani okolí.

Obecné nebezpečí	4	Nebezpečí zásahu el. proudem
Biologické nebezpečí		Nebezpečí pořezání
Nebezpečí v souvislosti s hořlavými materiály	i	Upozorňuje na důležité informace, které se nevztahují k nebezpečím.
Nebezpečí zranění prstů nebo rukou při neúmyslném zavření mechanických dílů.		

Bezpečnostní pokyny



Při nedodržení těchto bezpečnostních pokynů mohou vznikat nebezpečné situace, které mohou vést k těžkým nebo smrtelným zraněním osob, jestliže nezabráníte jejich vzniku.

- VÝSTRAHA
- Dodržujte bezpečnostní pokyny.
- Míchadlo je dovoleno používat jen v souladu s jeho určením. Použití v rozporu s určením může vést ke škodám, kontaminaci a zraněním se smrtelnými následky.
- Míchadlo smí provozovat jen proškolený personál.
- Provozovatel je povinen zajistit používání vhodných osobních ochranných prostředků. Dodržujte doporučení Příručky o biologické bezpečnosti v laboratořích (Laboratory Biosafety Manual zpracovaný Světovou zdravotnickou organizací WHO) a další platná národní doporučení.



Škody v důsledku nesprávného elektrického napájení.

Zkontrolujte, že je míchadlo připojeno jen do řádně uzemněné elektrické zásuvky.

VÝSTRAHA



Nebezpečí při zacházení s nebezpečnými látkami.

VÝSTRAHA

- Při práci s korozivními vzorky (roztoky solí, kyseliny, zásady) pečlivě vyčistěte míchadlo i díly příslušenství.
- Míchadlo není inertizované ani chráněné proti výbuchu. Míchadlo nikdy nepoužívejte ve výbušném prostředí.
- Nemíchejte toxické materiály, radioaktivní materiály ani patogenní mikroorganizmy bez provedení vhodných bezpečnostních opatření.
- Při míchání nebezpečných materiálů dodržujte doporučení Příručky o biologické bezpečnosti v laboratořích (Laboratory Biosafety Manual zpracovaný Světovou zdravotnickou organizací WHO) a další platná národní doporučení. Při míchání mikrobiologických vzorků rizikové skupiny II (dle Příručky o biologické bezpečnosti v laboratořích zpracované Světovou zdravotnickou organizací WHO) je nutné používat aerosoltěsná biotěsnění. Příručku o biologické bezpečnosti v laboratořích zpracované Světovou zdravotnickou organizací WHO) je nutné používat aerosoltěsná biotěsnění. Příručku o biologické bezpečnosti v laboratořích Laboratory Biosafety Manual najdete na internetové stránce Světové zdravotnické organizace (www.who.int). U materiálů vyšší rizikové skupiny je nutné provést další bezpečnostní opatření.
- Po kontaminaci míchadla nebo dílů toxiny nebo patogenními látkami je nutné provést vhodná dezinfekční a dekontaminační opatření ("Dekontaminace" na straně 124, "Dezinfekce" na straně 124).
- Nastane-li nebezpečná situace, přerušte napájení míchadla a ihned opusťte prostor.



VÝSTRAHA

Poškození zdraví infekčními látkami.

Proniknou-li nedopatřením kapaliny nebo jiné materiály pod základnu, ihned míchadlo vypněte, vytáhněte zástrčku z elektrické sítě a sejměte základnu ("Základny" na straně 18).

Utřete rozlitou kapalinu v souladu se standardním laboratorním postupem. Použijte vhodné osobní ochranné prostředky.



VÝSTRAHA

Poškození zdraví při míchání výbušných nebo hořlavých materiálů či látek.

Nemíchejte výbušné ani hořlavé materiály či látky.



Popálení na horkém povrchu.

POZOR

V míchadle s regulací teploty se mohou povrchy (např. základna a díly příslušenství) pod krytem velmi zahřívat. Nedotýkejte se horkých povrchů. Počkejte, než horké povrchy vychladnou.



POZOR

Nebezpečí pořezání skleněnými střepy.

Uvolněná základna a části příslušenství, např. držáky baněk, mohou při vypadnutí nádob z míchadla způsobit prasknutí skla.

Zkontrolujte, že jsou základna i části příslušenství řádně nainstalované, použijte správné nářadí a šrouby. Zkontrolujte, že jsou držáky baněk při zvolených otáčkách s to udržet nádoby/vzorky dané hmotnosti. Držáky baněk s individuálním nákladem mohou být případně vhodné jen pro otáčky ležící pod mezními hodnotami stability míchadla. Další informace k tomuto tématu najdete v oddíle "3. 4. Nakládání a použití v souladu s určením" na straně 115.

Věnujte pozornost neobvyklým zvukům. Ty mohou indikovat uvolněnou základnu nebo uvolněné části příslušenství.



Poškození míchadla nebo chybná funkce v důsledku poškození ovládacího panelu.

- Neprovozujte míchadlo s poškozeným ovládacím panelem.
- Vypněte míchadlo. Vytáhněte zástrčku z elektrické sítě. Nechte ovládací panel vyměnit autorizovaným servisním technikem.



POZOR

Biologické ohrožení rozbitými nebo netěsnými nádobami.

Neodborně nainstalované příslušenství může vést k rozlití vzorků.

- Zkontrolujte, že jsou díly příslušenství řádně nainstalované, použijte správné nářadí a šrouby.
- Zkontrolujte, že díly příslušenství řádně sedí na základně.
- Nádoby používejte vždy jen s příslušenstvím, které odpovídá jejich velikosti.
- Nádoby je nutné instalovat řádně a v neporušeném stavu.



POZOR

Neodborné naložení a poškozené příslušenství mohou ohrozit bezpečnost.

- Vždy dbejte na maximálně rovnoměrné rozložení nákladu (příslušenství a vzorků), především při provozu dvoupatrové základny.
- Nepoužívejte díly příslušenství vykazující známky koroze nebo praskliny. Pro další informace kontaktujte zákaznický servis.
- Používejte jen správně naložené míchadlo.
- Míchadlo nikdy nepřetěžujte.
- Před uvedením míchadla do provozu zkontrolujte, že je příslušenství řádně nainstalované. Postupujte podle pokynů v oddíle "Příslušenství" na straně 88.

POZOR



POZOR

Nebezpečí zranění při nedodržení provozních zásad.

- Míchadlo neprovozujte bez řádně namontované základny.
- Míchadlo nepoužívejte s poškozenými nebo demontovanými díly pláště.
- Míchadlo nikdy neprovozujte s vadnou pružinou krytu. Vadná pružina krytu není s to udržet kryt bezpečně ve zcela otevřené pozici.
- S míchadlem v provozu nehýbejte.
- O míchadlo se neopírejte.
- Míchadlo nakládejte a vykládejte vždy až po jeho úplném zastavení, úplné zastavení je indikováno na ovládacím panelu.
- Na míchadlo v provozu nic nepokládejte.
- Nedotýkejte se základny ani částí příslušenství, jestliže je míchadlo v provozu.
- Personál obsluhy nesmí otevírat plášť míchadla.



Nekompatibilní části příslušenství mohou ohrozit bezpečnost.

Používejte v tomto míchadle jen příslušenství schválené společností Thermo Fisher Scientific. Aktuální seznamy najdete na <u>www.thermofisher.</u> <u>com</u>.

UPOZORNĚNÍ



Vypnutí míchadla:

Stiskněte tlačítko STOP. Vypněte míchadlo hlavním vypínačem. Vytáhněte zástrčku z elektrické sítě. V případě nouze přerušte přívod proudu.

UPOZORNĚNÍ

1. Technické specifikace

1.1. Technické údaje

Míchadlo Thermo Scientific Solaris 2000				
Rozsah otáček	15–525 ot./min			
Procesní čas	99 h 59 m (v krocích po 1 minutě) nebo trvalý provoz			
Hladina hluku při max. otáčkách	50 dB (A) (1 m před zařizením ve výšce 1,6 m)			
Max. náklad (vč. základny, příslušenství a vzorků)	25 kg (55 lb)			
Elektrické připojení	100–240 V ±10 %, 50 / 60 Hz			
Příkon	60 W			
Podmínky okolí				
Skladování a přeprava	Teplota: -10 °C až 55 °C Vlhkost: 15 až 85 %			
Při provozu	Použití ve vnitřních prostorech Výška do 3000 m nad mořem Teplota: 5 až 40 °C Max. relativní vlhkost 80 % do 31 °C, lineární pokles až na 50 % rel. vlhkost při 40 °C.			
Stupeň znečištění	2			
Přepěťová kategorie	Ш			
lb	20			
Rozhraní	USB	Ethernet		
	2x USB-A 2.0	RJ45		
Rozměry	Délka	Šířka	Výška	
	47 cm (18,5 palců)	37 cm (14,5 palců)	15 cm (5,5 palců)	
Hmotnost	20,9 kg (46,0 lb)			

Tabulka 1: Technické údaje Solaris 2000

Technické údaje

Míchadlo Thermo Scientific Solaris 4000

Rozsah otáček	15–525 ot./min			
Procesní čas	99 h 59 m (v krocích po 1 minutě) nebo trvalý provoz			
Hladina hluku při max. otáčkách	54 dB (A) (1 m před zařízením ve výšce 1,6 m)			
Max. náklad (vč. základny, příslušenství a vzorků)	43 kg (95 lb)			
Elektrické připojení	100–240 V ±10 %, 50 / 60 Hz			
Příkon	80 W			
Podmínky okolí				
Skladování a přeprava	Teplota: -10 °C až 55 °C Vlhkost: 15 až 85 %			
Při provozu	Použití ve vnitřních prostore Výška do 3000 m nad moře Teplota: 5 až 40 °C Max. relativní vlhkost 80 % d	ch m o 31 °C, lineární pokles až na :	50 % rel. vlhkost při 40 °C.	
Stupeň znečištění	2			
Přepěťová kategorie	11			
IP	20			
Rozhraní	USB	Ethernet		
	2x USB-A 2.0	RJ45		
Rozměry	Délka	Šířka	Výška	
	65 cm (25,6 palců)	58 cm (22,8 palců)	18 cm (7 palců)	
Hmotnost	75,1 kg (165,5 lb)			

Tabulka 2: Technické údaje Solaris 4000

Míchadlo Thermo Scientific Solaris 2000 I

Rozsah otáček	15–525 ot./min			
Nastavovací rozsah teploty	30-60 °C			
Teplotní stabilita v baňce při 37 °C	Teplota okolí 23 °C; za 1 hod. při stabilní teplotě v komoře \pm 0,1 °C			
Teplotní rovnoměrnost v baňce při 37 °C	Teplota okolí 23 °C; za 1 hod. při stabilní teplotě v komoře \pm 0,5 °C			
Procesní čas	99 h 59 m (v krocích po 1 minutě) nebo trvalý provoz			
Hladina hluku při max. otáčkách	52 dB (A) (1 m před zařízením ve výšce 1,6 m)			
Max. náklad (vč. základny, příslušenství a vzorků)	16 kg (35 lb), vč. základny 11x14 se 2,3 kg (5,1 lb)			
Elektrické připojení	100–120, 200–240 V ±10 %	, 50 / 60 Hz		
Příkon	900 W			
Podmínky okolí				
Skladování a přeprava	Teplota: -10 °C až 55 °C Vlhkost: 15 až 85 %			
Při provozu	Použití ve vnitřních prostorech Výška do 3000 m nad mořem Teplota: 5 až 40 °C Max. relativní vlhkost 80 % do 31 °C, lineární pokles až na 50 % rel. vlhkost při 40 °C.			
	Vyska do 3000 m nad morer Teplota: 5 až 40 °C Max. relativní vlhkost 80 % do	n o 31 °C, lineární pokles až na {	50 % rel. vlhkost při 40 °C.	
Stupeň znečištění	Vyska do 3000 m nad morer Teplota: 5 až 40 °C Max. relativní vlhkost 80 % do 2	n o 31 °C, lineární pokles až na š	50 % rel. vlhkost při 40 °C.	
Stupeň znečištění Přepěťová kategorie	Vyska do 3000 m nad morer Teplota: 5 až 40 °C Max. relativní vlhkost 80 % do 2 II	n o 31 °C, lineární pokles až na t	50 % rel. vlhkost při 40 °C.	
Stupeň znečištění Přepěťová kategorie IP	Vyska do 3000 m nad morer Teplota: 5 až 40 °C Max. relativní vlhkost 80 % do 2 II 20	n o 31 °C, lineární pokles až na t	50 % rel. vlhkost při 40 °C.	
Stupeň znečištění Přepěťová kategorie IP Rozhraní	Vyska do 3000 m had morer Teplota: 5 až 40 °C Max. relativní vlhkost 80 % do 2 II 20 USB	n o 31 °C, lineární pokles až na t Ethernet	50 % rel. vihkost při 40 °C.	
Stupeň znečištění Přepěťová kategorie IP Rozhraní	Vyska do 3000 m had morer Teplota: 5 až 40 °C Max. relativní vlhkost 80 % d 2 II 20 USB 2x USB-A 2.0	n o 31 °C, lineární pokles až na 5 Ethernet RJ45	50 % rel. vlhkost při 40 °C.	
Stupeň znečištění Přepěťová kategorie IP Rozhraní Rozměry	Vyska do 3000 m nad morer Teplota: 5 až 40 °C Max. relativní vlhkost 80 % d 2 II 20 USB 2x USB-A 2.0 Délka	n o 31 °C, lineární pokles až na 5 Ethernet RJ45 Šířka	50 % rel. vlhkost při 40 °C. Výška	
Stupeň znečištění Přepěťová kategorie IP Rozhraní Rozměry	Vyska do 3000 m had morer Teplota: 5 až 40 °C Max. relativni vlhkost 80 % do 2 II 20 USB 2x USB-A 2.0 Délka 70 cm (27,6 palců)	n o 31 °C, lineární pokles až na 8 Ethernet RJ45 Šiřka 36 cm (14,2 palců)	50 % rel. vlhkost při 40 °C. Výška 46 cm (18,1 palců)	
Stupeň znečištění Přepěťová kategorie IP Rozhraní Rozměry S otevřeným krytem	Vyska do 3000 m nad morer Teplota: 5 až 40 °C Max. relativní vlhkost 80 % d 2 II 20 USB 2x USB-A 2.0 Délka 70 cm (27,6 palců)	n o 31 °C, lineární pokles až na 5 Ethernet RJ45 Šířka 36 cm (14,2 palců)	50 % rel. vlhkost při 40 °C. Výška 46 cm (18,1 palců) 79 cm (30,8 palců)	

Tabulka 3: Technické údaje Solaris 2000 l

Míchadlo Thermo Scientific Solaris 4000 I



Tabulka 4: Technické údaje Solaris 4000 I

Míchadlo Thermo Scientific Solaris 2000 R

Rozsah otáček	15–525 ot./min			
Nastavovací rozsah teploty	5–60 °C			
Teplotní stabilita v baňce při 37 °C	Teplota okolí 23 °C; za 1 hod. při stabilní teplotě v komoře \pm 0,1 °C			
Teplotní rovnoměrnost v baňce při 37 °C	Teplota okolí 23 °C; za 1 hod. při stabilní teplotě v kornoře \pm 0,5 °C			
Procesní čas	99 h 59 m (v krocích po 1 minutě) nebo trvalý provoz			
Hladina hluku při max. otáčkách	56 dB (A) (1 m před zařízen	ím ve výšce 1,6 m)		
Max. náklad (vč. základny, příslušenství a vzorků)	16 kg (35 lb), vč. základny 11x14 se 2,3 kg (5,1 lb)			
Elektrické připojení	100-240 V ±10 %, 50 / 60 H	łz		
Příkon	350 W			
Podmínky okolí				
Skladování a přeprava	Teplota: -10 °C až 55 °C Vlhkost: 15 až 85 %			
Při provozu	Použití ve vnitřních prostorech Výška do 3000 m nad mořem Teplota: 5 až 40 °C Max. relativní vlhkost 80 % do 31 °C, lineární pokles až na 50 % rel. vlhkost při 40 °C.			
Stupeň znečištění	2			
Stupeň znečištění Přepěťová kategorie	2			
Stupeň znečištění Přepěťová kategorie IP	2 II 20			
Stupeň znečištění Přepěťová kategorie IP Rozhraní	2 II 20 USB	Ethernet		
Stupeň znečištění Přepěťová kategorie IP Rozhraní	2 II 20 USB 2x USB-A 2.0	Ethernet RJ45		
Stupeň znečištění Přepěťová kategorie IP Rozhraní Rozměry	2 II 20 USB 2x USB-A 2.0 Délka	Ethernet RJ45 Šířka	Výška	
Stupeň znečištění Přepěťová kategorie IP Rozhraní Rozměry	2 II 20 USB 2x USB-A 2.0 Délka 70 cm (27,6 palců)	Ethernet RJ45 Šiřka 36 cm (14,2 palců)	Výška 46 cm (18,1 palců)	
Stupeň znečištění Přepěťová kategorie IP Rozhraní Rozměry S otevřeným krytem	2 II 20 USB 2x USB-A 2.0 Délka 70 cm (27,6 palců)	Ethernet RJ45 Šiřka 36 cm (14,2 palců)	Výška 46 cm (18,1 palců) 79 cm (30,8 palců)	

Tabulka 5: Technické údaje Solaris 2000 R

Míchadlo Thermo Scientific Solaris 4000 R

Rozsah otáček	15–525 ot./min		
Nastavovací rozsah teploty	4–60 °C		
Teplotní stabilita v baňce při 37 °C	Teplota okolí 23 °C; za 1 hod. při stabilní teplotě v komoře \pm 0,1 °C		
Teplotní rovnoměrnost v baňce při 37 °C	Teplota okolí 23 °C; za 1 hod. při stabilní teplotě v komoře \pm 0,5 °C		
Procesní čas	99 h 59 m (v krocích po 1 minutě) nebo trvalý provoz		
Hladina hluku při max. otáčkách	56 dB (A) (1 m před zařízenír	n ve výšce 1,6 m)	
Max. náklad (vč. základny, příslušenství a vzorků)	32 kg (70,5 lb), vč. základny 18x18 se 4,3 kg (9,5 lb)		
Elektrické připojení	100–240 V ±10 %, 50 / 60 H;	z	
Příkon	750 W		
Podmínky okolí			
Skladování a přeprava	Teplota: -10 °C až 55 °C Vlhkost: 15 až 85 %		
Při provozu	Použití ve vnitřních prostorec Výška do 3000 m nad mořem Teplota: 5 až 40 °C Max. relativní vlhkost 80 % do	h o 31 °C, lineární pokles až na 5	50 % rel. vlhkost při 40 °C.
Stupeň znečištění	2		
Přepěťová kategorie	II		
IP	20		
Rozhraní	USB	Ethernet	
	2x USB-A 2.0	RJ45	
Rozměry	Délka	Šířka	Výška
	77 cm (30,3 palců)	57 cm (22,4 palců)	55 cm (21,7 palců)
S otevřeným krytem			94 cm (36,7 palců)
Hmotnost	74,9 kg (165,1 lb)		

Tabulka 6: Technické údaje Solaris 4000 R

Příslušenství

1.2. Příslušenství



Nekompatibilní části příslušenství mohou ohrozit bezpečnost.

Používejte v tomto míchadle jen příslušenství schválené společností Thermo Fisher Scientific.

POZOR

Aktuální seznamy najdete na www.thermofisher.com.

Příslušenství

1. 2. 1. Základny

Základna		
Solaris 2000	Hmotnost	7
Univerzální základna Thermo Scientific Solaris 12x14	2,7 kg (5,9 lb)	SK1214
Dvoupatrová univerzální základna Thermo Scientific Solaris 12x14	6,4 kg (14,1 lb)	SK1214D
Sada pro upgrade, dvoupatrová univerzální základna Thermo Scientific Solaris 12x14		SK1214DK
Univerzální základna Thermo Scientific Solaris 18x18	4,3 kg (9,4 lb)	SK1818
Dvoupatrová univerzální základna Thermo Scientific Solaris 18x18	9,7 kg (21,3 lb)	SK1818D
Sada pro upgrade, dvoupatrová univerzální základna Thermo Scientific Solaris 18x18		SK1818DK
Univerzální základna Thermo Scientific Solaris 18x24	6,0 kg (13,2 lb)	SK1824
Solaris 4000		
Univerzální základna Thermo Scientific Solaris 18x30	7,5 kg (16,5 lb)	SK1830
Dvoupatrová univerzální základna Thermo Scientific Solaris 18x30	16,3 kg (35,9 lb)	SK1830D
Sada pro upgrade, dvoupatrová univerzální základna Thermo Scientific Solaris 18x30		SK1830DK
Univerzální základna Thermo Scientific Solaris 36x24	12,0 kg (26,4 lb)	SK3624
Solaris 2000 I / 2000 R		
Univerzální základna Thermo Scientific Solaris 11x14	2,3 kg (5,1 lb)	SK1114
Solaris 4000 I / 4000 R		
Univerzální základna Thermo Scientific Solaris 18x18	4,3 kg (9,5 lb)	SK1818
Doplňující sady a příslušenství		
Sada náhradních dílů pro držák baněk (šrouby)		SK1001
Sada náhradních dílů pro základnu, malá (šrouby základny, nářadí)		SK0100
Sada náhradních dílů pro základnu, velká (šrouby základny, nářadí)		SK0101
Śroubovák pro dily příslušenství		75004131

Tabulka 7: Nabízené základny

1. 2. 2. Držáky baněk



Pro základny Solaris 2000

	Č. výrobku	Univerzální			Dvoup	atrová
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Erlenmeyerova baňka 10 ml	30150BI	72	113	157	140	226
Erlenmeyerova baňka 25 ml	30151	42	64	80	80	124
Erlenmeyerova baňka 50 ml	30152BI	42	64	80	80	124
Erlenmeyerova baňka 125 ml	30153	15	32	40	30	52
Erlenmeyerova baňka 250 ml	30154BI	9	16	24	16	32
Erlenmeyerova baňka 300 ml	30155	9	16	20	16	32
Erlenmeyerova baňka 500 ml	30156BI	9	16	20	16	32
Erlenmeyerova baňka 1 l	30157BI	4	9	10	8	16
Erlenmeyerova baňka 2 l	30158	3	5	6	-	-
Erlenmeyerova baňka 4 l	30159	1	4	4	-	-
Erlenmeyerova baňka 5 l	30159B	1	2	4	-	-
Erlenmeyerova baňka 6 l	30160	1	2	2	-	-
Fernbachova baňka 2800 ml	30162	1	4	4	-	-
Kultivační baňka 2500 ml, nízká	30161	1	1	2	-	-

Tabulka 8: Nabízené držáky baněk pro základny Solaris 2000

Pro základny Solaris 4000

	Č. výrobku	Univerzální		Dvoupatrová
		18 x 30	36 x 24	18 x 30
Erlenmeyerova baňka 10 ml	30150BI	203	187	402
Erlenmeyerova baňka 25 ml	30151	112	187	220
Erlenmeyerova baňka 50 ml	30152BI	112	187	220
Erlenmeyerova baňka 125 ml	30153	46	83	92
Erlenmeyerova baňka 250 ml	30154BI	28	40	56
Erlenmeyerova baňka 300 ml	30155	28	40	56
Erlenmeyerova baňka 500 ml	30156BI	28	40	56
Erlenmeyerova baňka 1 l	30157BI	14	20	28
Erlenmeyerova baňka 2 l	30158	6	11	12
Erlenmeyerova baňka 4 l	30159	6	8	-
Erlenmeyerova baňka 5 l	30159B	5	8	-
Erlenmeyerova baňka 6 l	30160	3	6	-
Fernbachova baňka 2800 ml	30162	6	8	12
Kultivační baňka 2500 ml, nízká	30161	3	6	6

Tabulka 9: Nabízené držáky baněk pro základny Solaris 4000

Pro základny Solaris 2000 I / 2000 R

	Č. výrobku	Univerzální
		11 x 14
Erlenmeyerova baňka 10 ml	30150BI	59
Erlenmeyerova baňka 25 ml	30151	35
Erlenmeyerova baňka 50 ml	30152BI	35
Erlenmeyerova baňka 125 ml	30153	15
Erlenmeyerova baňka 250 ml	30154BI	7
Erlenmeyerova baňka 300 ml	30155	7
Erlenmeyerova baňka 500 ml	30156BI	7
Erlenmeyerova baňka 1 l	30157BI	4
Erlenmeyerova baňka 2 l	30158	2
Erlenmeyerova baňka 4 l	30159	-
Erlenmeyerova baňka 5 l	30159B	-
Erlenmeyerova baňka 6 l	30160	-
Fernbachova baňka 2800 ml	30162	1
Kultivační baňka 2500 ml, nízká	30161	1

Tabulka 10: Nabízené držáky baněk pro základny Solaris 2000 I / 2000 R

Pro základny Solaris 4000 I / 4000 R

	Č. výrobku	Univerzální
		18 x 18
Erlenmeyerova baňka 10 ml	30150BI	113
Erlenmeyerova baňka 25 ml	30151	64
Erlenmeyerova baňka 50 ml	30152BI	64
Erlenmeyerova baňka 125 ml	30153	32
Erlenmeyerova baňka 250 ml	30154BI	16
Erlenmeyerova baňka 300 ml	30155	16
Erlenmeyerova baňka 500 ml	30156BI	16
Erlenmeyerova baňka 1 l	30157BI	9
Erlenmeyerova baňka 2 l	30158	6
Erlenmeyerova baňka 4 l	30159	-
Erlenmeyerova baňka 5 l	30159B	-
Erlenmeyerova baňka 6 l	30160	-
Fernbachova baňka 2800 ml	30162	4
Kultivační baňka 2500 ml, nízká	30161	1

Tabulka 11: Nabízené držáky baněk pro základny Solaris 4000 I / 4000 R

1. 2. 3. Držáky baněk pro mikrotitrační/DeepWell destičky



Pro základny Solaris 2000

	Č. výrobku	Univerzální		Dvoup	atrová	
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Mikrotitrační/DeepWell destička	30175	7	12	16	14	24

Tabulka 12: Kapacita držáků pro mikrotitrační/DeepWell destičky pro základny Solaris 2000

Pro základny Solaris 2000 I / 2000 R

	Č. výrobku	Univerzální
		11 x 14
Mikrotitrační/DeepWell destička	30175	6

Tabulka 13: Kapacita držáků pro mikrotitrační/DeepWell destičky pro základny Solaris 2000 I / 2000 R

Pro základny Solaris 4000

	Č. výrobku	Univerzální		Dvoupatrová
		18 x 30	36 x 24	18 x 30
Mikrotitrační/DeepWell destička	30175	22	36	42

Tabulka 14: Kapacita držáků pro mikrotitrační/DeepWell destičky pro základny Solaris 4000

Pro základny Solaris 4000 I / 4000 R

	Č. výrobku	Univerzální
		18 x 18
Mikrotitrační/DeepWell destička	30175	10

Tabulka 15: Kapacita držáků pro mikrotitrační/DeepWell destičky pro základny Solaris 4000 I / 4000 R

1. 2. 4. Stojany na zkumavky

Pro základny Solaris 2000



	Č. výrobku	Univerzální			Dvoupatrová	
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Poloviční velikost						
10–13 mm, červený, pole 6x6	30181	8	12	15	14	24
14–16 mm, oranžový, pole 6x6	30183	5	9	11	9	17
17–20 mm, bílý, pole 4x5	30185	7	11	14	13	21
21–25 mm, modrý, pole 4x4	30187	6	9	11	10	17
26–30 mm, zelený, pole 3x3	30189	6	9	12	11	18
pro mikrotitrační zkumavky, 1,5 ml, modrý, pole 4x6	30191	6	10	13	11	19
Plná velikost						
10-13 mm, pole 6x12	30180BI	3	7	10	6	13
14-16 mm, pole 6x12	30182	3	4	6	6	8
17–20 mm, pole 4x10	30184	3	5	7	6	8
21–25 mm, pole 4x10	30186	2	3	5	4	6
26–30 mm, pole 3x8	30188	3	4	6	5	8
pro mikrotitrační zkumavky, 1,5 ml, pole 8x12	30190	3	5	7	6	9

Tabulka 16: Nabízené stojany na zkumavky pro základny Solaris 2000

Příslušenství

Pro základny Solaris 4000

	Č. výrobku	Univerzální		Dvoupatrová
		18 x 30	36 x 24	18 x 30
Poloviční velikost				
10–13 mm, červený, pole 6x6	30181	21	32	40
14–16 mm, oranžový, pole 6x6	30183	14	20	26
17–20 mm, bílý, pole 4x5	30185	18	20	34
21–25 mm, modrý, pole 4x4	30187	14	22	25
26–30 mm, zelený, pole 3x3	30189	15	24	30
pro mikrotitrační zkumavky, 1,5 ml, modrý, pole 4x6	30191	18	24	34
Plná velikost				
10-13 mm, pole 6x12	30180BI	14	20	26
14-16 mm, pole 6x12	30182	9	12	16
17-20 mm, pole 4x10	30184	9	15	18
21-25 mm, pole 4x10	30186	7	9	13
26-30 mm, pole 3x8	30188	7	10	13
pro mikrotitrační zkumavky, 1,5 ml, pole 8x12	30190	9	12	17

Tabulka 17: Nabízené stojany na zkumavky pro základny Solaris 4000

Pro základny Solaris 2000 I / 2000 R

	Č. výrobku	Univerzální
		11 x 14
Poloviční velikost		
10–13 mm, červený, pole 6x6	30181	6
14–16 mm, oranžový, pole 6x6	30183	2
17–20 mm, bílý, pole 4x5	30185	4
21–25 mm, modrý, pole 4x4	30187	2
26–30 mm, zelený, pole 3x3	30189	4
pro mikrotitrační zkumavky, 1,5 ml, modrý, pole 4x6	30191	3
Plná velikost		
10-13 mm, pole 6x12	30180BI	3
14–16 mm, pole 6x12	30182	2
17–20 mm, pole 4x10	30184	2
21–25 mm, pole 4x10	30186	1
26–30 mm, pole 3x8	30188	2
pro mikrotitrační zkumavky, 1,5 ml, pole 8x12	30190	2

Tabulka 18: Nabízené stojany na zkumavky pro základny Solaris 2000 I / 2000 R

Pro základny Solaris 4000 I / 4000 R

	Č. výrobku	Univerzální
		18 x 18
Poloviční velikost		
10–13 mm, červený, pole 6x6	30181	12
14–16 mm, oranžový, pole 6x6	30183	8
17–20 mm, bílý, pole 4x5	30185	10
21–25 mm, modrý, pole 4x4	30187	8
26–30 mm, zelený, pole 3x3	30189	9
pro mikrotitrační zkumavky, 1,5 ml, modrý, pole 4x6	30191	8
Plná velikost		
10-13 mm, pole 6x12	30180BI	7
14-16 mm, pole 6x12	30182	4
17-20 mm, pole 4x10	30184	5
21-25 mm, pole 4x10	30186	3
26-30 mm, pole 3x8	30188	4
pro mikrotitrační zkumavky, 1,5 ml, pole 8x12	30190	5

Tabulka 19: Nabízené stojany na zkumavky pro základny Solaris 4000 I / 4000 R

1. 2. 5. Držáky stojanů na zkumavky s nastavitelným úhlem



Pro základny Solaris 2000

	Č. výrobku		Univerzáln	í	Dvoup	atrová
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
10-13 mm, pro 72 zkumavek	236090	3	4	6	6	8
16-20 mm, pro 40 zkumavek	236091	3	4	6	5	8
21-25 mm, pro 40 zkumavek	236092	3	3	4	5	6
26-30 mm, pro 24 zkumavek	236093	3	3	4	5	6

Tabulka 20: Nabízené držáky stojanů na zkumavky s nastavitelným úhlem pro základny Solaris 2000

Pro základny Solaris 4000

	Č. výrobku	Univerzální		Dvoupatrová
		18 x 30	36 x 24	18 x 30
10-13 mm, pro 72 zkumavek	236090	7	14	14
16-20 mm, pro 40 zkumavek	236091	7	14	14
21-25 mm, pro 40 zkumavek	236092	5	9	10
26-30 mm, pro 24 zkumavek	236093	6	9	12

Tabulka 21: Nabízené držáky stojanů na zkumavky s nastavitelným úhlem pro základny Solaris 4000

Pro základny Solaris 2000 I / 2000 R

	Č. výrobku	Univerzální
		11 x 14
10-13 mm, pro 72 zkumavek	236090	2
16-20 mm, pro 40 zkumavek	236091	2
21-25 mm, pro 40 zkumavek	236092	1
26-30 mm, pro 24 zkumavek	236093	1

Tabulka 22: Nabízené držáky stojanů na zkumavky s nastavitelným úhlem pro základny Solaris 2000 l / 2000 R

Pro základny Solaris 4000 I / 4000 R

	Č. výrobku	Univerzální
		18 x 18
10-13 mm, pro 72 zkumavek	236090	4
16-20 mm, pro 40 zkumavek	236091	3
21-25 mm, pro 40 zkumavek	236092	3
26-30 mm, pro 24 zkumavek	236093	3

Tabulka 23: Nabízené držáky stojanů na zkumavky s nastavitelným úhlem pro základny Solaris 4000 I / 4000 R

1. 2. 6. Nastavitelné držáky nádob



Pro základny Solaris 2000

	Č. výrobku	I	Univerzáln	í	Dvoup	atrová
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Jednostupňový výškově nastavitelný držák nádob 11"x14"	75004104	-	-	-	-	-
Dvoustupňový výškově nastavitelný držák nádob 12"x14"	75004102	1	1	1	-	2
Dvoustupňový výškově nastavitelný držák nádob 18"x18"	75004103	-	1	1	-	-
Nastavitelný držák jedné nádoby	75004101	4	4	6	6	8

Tabulka 24: Nabízené nastavitelné držáky nádob pro základny Solaris 2000

Pro základny Solaris 4000

	Č. výrobku	Univerzální		Dvoupatrová
		18 x 30	36 x 24	18 x 30
Jednostupňový výškově nastavitelný držák nádob 11"x14"	75004104	-	-	-
Dvoustupňový výškově nastavitelný držák nádob 12"x14"	75004102	2	2	4
Dvoustupňový výškově nastavitelný držák nádob 18"x18"	75004103	1	2	2
Nastavitelný držák jedné nádoby	75004101	8	12	16

Tabulka 25: Nastavitelné držáky nádob pro základny Solaris 4000

Pro základny Solaris 2000 I / 2000 R

	Č. výrobku	Univerzální
		11 x 14
Jednostupňový výškově nastavitelný držák nádob 11"x14"	75004104	1
Dvoustupňový výškově nastavitelný držák nádob 12"x14"	75004102	-
Dvoustupňový výškově nastavitelný držák nádob 18"x18"	75004103	-
Nastavitelný držák jedné nádoby	75004101	2

Tabulka 26: Nastavitelné držáky nádob pro základny Solaris 2000 I / 2000 R

Pro základny Solaris 4000 I / 4000 R

	Č. výrobku	Univerzální
		18 x 18
Jednostupňový výškově nastavitelný držák nádob 11"x14"	75004104	-
Dvoustupňový výškově nastavitelný držák nádob 12"x14"	75004102	1
Dvoustupňový výškově nastavitelný držák nádob 18"x18"	75004103	1
Nastavitelný držák jedné nádoby	75004101	4

Tabulka 27: Nastavitelné držáky nádob pro základny Solaris 4000 I / 4000 R

1. 2. 7. Čtvercové držáky nádob



Pro základny Solaris 2000

	Č. výrobku	I	Univerzáln	í	Dvoup	atrová
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Čtvercový držák lahve, 125 ml	75004106	12	16	24	21	32
Čtvercový držák lahve, 250 ml	75004107	9	16	22	18	32
Čtvercový držák lahve, 500 ml	75004108	6	9	12	11	18
Čtvercový držák lahve, 1000 ml	75004109	4	9	11	8	18

Tabulka 28: Nabízené čtvercové držáky nádob pro základny Solaris 2000

Pro základny Solaris 4000

	Č. výrobku	Univerzální		Dvoupatrová
		18 x 30	36 x 24	18 x 30
Čtvercový držák lahve, 125 ml	75004106	28	54	56
Čtvercový držák lahve, 250 ml	75004107	28	52	56
Čtvercový držák lahve, 500 ml	75004108	15	24	30
Čtvercový držák lahve, 1000 ml	75004109	15	24	28

Tabulka 29: Nabízené čtvercové držáky nádob pro základny Solaris 4000

Pro základny Solaris 2000 I / 2000 R

	Č. výrobku	Univerzální
		11 x 14
Čtvercový držák lahve, 125 ml	75004106	9
Čtvercový držák lahve, 250 ml	75004107	7
Čtvercový držák lahve, 500 ml	75004108	3
Čtvercový držák lahve, 1000 ml	75004109	2

Tabulka 30: Nabízené čtvercové držáky nádob pro základny Solaris 2000 I / 2000 R

Pro základny Solaris 4000 I / 4000 R

	Č. výrobku	Univerzální	
		18 x 18	
Čtvercový držák lahve, 125 ml	75004106	16	
Čtvercový držák lahve, 250 ml	75004107	16	
Čtvercový držák lahve, 500 ml	75004108	9	
Čtvercový držák lahve, 1000 ml	75004109	9	

Tabulka 31: Nabízené čtvercové držáky nádob pro základny Solaris 4000 I / 4000 R

1. 2. 8. Stojany pro kádinky Nalgene

Pro základny Solaris 2000



	Č. výrobku	Univerzální		Dvoup	Dvoupatrová	
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Velikost 11x14 / 12x14			S	tojany / místa		
30 ml kádinka Nalgene, 36 míst	75004116	1 / 36	-	-	2/72	-
50 ml kádinka Nalgene, 28 míst	75004129	1 / 28	-	-	2 / 56	-
100 ml kádinka Nalgene, 16 míst	75004118	1 / 16	-	-	2/32	-
250 ml kádinka Nalgene, 9 míst	75004119	1/9	-	-	2 / 18	-
400 ml kádinka Nalgene, 8 míst	75004120	1/8	-	-	2/16	-
600 ml kádinka Nalgene, 5 míst	75004121	1/5	-	-	2/10	-
Velikost 9x18			S	tojany / místa		
30 ml kádinka Nalgene, 32 míst	75004110	-	2 / 64	2/64	-	4 / 128
50 ml kádinka Nalgene, 24 míst	75004128	-	2 / 48	2 / 48	-	4 / 96
100 ml kádinka Nalgene, 15 míst	75004112	-	2 / 30	2/30	-	4 / 60
250 ml kádinka Nalgene, 8 míst	75004113	-	2 / 16	2 / 16	-	4 / 32
400 ml kádinka Nalgene, 6 míst	75004114	-	2/12	2 / 12	-	4 / 24
600 ml kádinka Nalgene, 4 míst	75004115	-	2/8	2/8	-	4 / 16

Tabulka 32: Nabízené stojany pro kádinky Nalgene pro základny Solaris 2000

Pro základny Solaris 4000

	Č. výrobku	Univerzální		Dvoupatrová
		18 x 30	36 x 24	18 x 30
Velikost 11x14 / 12x14			Stojany / místa	
30 ml kádinka Nalgene, 36 míst	75004116	-	-	-
50 ml kádinka Nalgene, 28 míst	75004129	-	-	-
100 ml kádinka Nalgene, 15 míst	75004118	-	-	-
250 ml kádinka Nalgene, 9 míst	75004119	-	-	-
400 ml kádinka Nalgene, 8 míst	75004120	-	-	-
600 ml kádinka Nalgene, 5 míst	75004121	-	-	-
Velikost 9x18			Stojany / místa	
30 ml kádinka Nalgene, 32 míst	75004110	3 / 96	4 / 128	6 / 192
50 ml kádinka Nalgene, 24 míst	75004128	3/72	4 / 96	6 / 144
100 ml kádinka Nalgene, 15 míst	75004112	3 / 45	4 / 60	6 / 90
250 ml kádinka Nalgene, 8 míst	75004113	3 / 24	4 / 32	6 / 48
400 ml kádinka Nalgene, 6 míst	75004114	3 / 18	4 / 24	6 / 36
600 ml kádinka Nalgene, 4 míst	75004115	3 / 12	4 / 16	6 / 24

Tabulka 33: Nabízené stojany pro kádinky Nalgene pro základny Solaris 4000

Pro základny Solaris 2000 I / 2000 R

	Č. výrobku	Univerzální
		11 x 14
Velikost 11x14 / 12x14		Stojany / místa
30 ml kádinka Nalgene, 36 míst	75004116	1/36
50 ml kádinka Nalgene, 28 míst	75004129	1 / 28
100 ml kádinka Nalgene, 16 míst	75004118	1 / 16
250 ml kádinka Nalgene, 9 míst	75004119	1/9
400 ml kádinka Nalgene, 8 míst	75004120	1/8
600 ml kádinka Nalgene, 5 míst	75004121	1/5
Velikost 9x18		Stojany / místa
30 ml kádinka Nalgene, 32 míst	75004110	-
50 ml kádinka Nalgene, 24 míst	75004128	-
100 ml kádinka Nalgene, 15 míst	75004112	-
250 ml kádinka Nalgene, 8 míst	75004113	-
400 ml kádinka Nalgene, 6 míst	75004114	-
600 ml kádinka Nalgene, 4 míst	75004115	-

Tabulka 34: Nabízené stojany pro kádinky Nalgene pro základny Solaris 2000 I / 2000 R
Pro základny Solaris 4000 I / 4000 R

	Č. výrobku	Univerzální
		18 x 18
Velikost 11x14 / 12x14		Stojany / místa
30 ml kádinka Nalgene, 36 míst	75004116	-
50 ml kádinka Nalgene, 28 míst	75004129	-
100 ml kádinka Nalgene, 16 míst	75004118	-
250 ml kádinka Nalgene, 9 míst	75004119	-
400 ml kádinka Nalgene, 8 míst	75004120	-
600 ml kádinka Nalgene, 5 míst	75004121	-
Velikost 9x18		Stojany / místa
30 ml kádinka Nalgene, 32 míst	75004110	2 / 64
50 ml kádinka Nalgene, 24 míst	75004128	2/48
100 ml kádinka Nalgene, 15 míst	75004112	2/30
250 ml kádinka Nalgene, 8 míst	75004113	2 / 16
400 ml kádinka Nalgene, 6 míst	75004114	2 / 12
600 ml kádinka Nalgene, 4 míst	75004115	2/8

Tabulka 35: Nabízené stojany pro kádinky Nalgene pro základny Solaris 4000 I / 4000 R

1. 2. 9. Držáky pro dělicí nálevky

Pro základny Solaris 2000

1. 2. 9. Dizaky più		y				10
Pro základny Solaris	<u>2000</u>				0	- Ho
	Č. výrobku		Univerzáln	í	Dvoup	atrová
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Jednobaňkový držák pro dělicí nálevku, 250 až 2000 ml	75004125	2	3	4	2	5
Držák pro dělicí nálevku, svislá montáž		4	4	6	-	-

Tabulka 36: Nabízené držáky pro dělicí nálevky pro základny Solaris 2000

Pro základny Solaris 4000

	Č. výrobku	Univerzální		Dvoupatrová
		18 x 30	36 x 24	18 x 30
Jednobaňkový držák pro dělicí nálevku, 250 až 2000 ml	75004125	5	7	9
Držák pro dělicí nálevku, svislá montáž		8	12	-

Tabulka 37: Nabízené držáky pro dělicí nálevky pro základny Solaris 4000

Pro základny Solaris 2000 I / 2000 R

	Č. výrobku	Univerzální
		11 x 14
Jednobaňkový držák pro dělicí nálevku, 250 až 2000 ml	75004125	-
Držák pro dělicí nálevku, svislá montáž		-

Tabulka 38: Nabízené držáky pro dělicí nálevky pro základny Solaris 2000 I / 2000 R

Pro základny Solaris 4000 I / 4000 R

	Č. výrobku	Univerzální	
		18 x 18	
Jednobaňkový držák pro dělicí nálevku, 250 až 2000 ml	75004125	-	
Držák pro dělicí nálevku, svislá montáž		-	

Tabulka 39: Nabízené držáky pro dělicí nálevky pro základny Solaris 4000 I / 4000 R

1. 2. 10. Univerzální odkládací tác

Pro základny Solaris 2000

	Č. výrobku	Univerzální			Dvoupatrová		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18	
Univerzální odkládací tác 11x14 / 12x14	75004123	1	-	-	2	-	
Univerzální odkládací tác 9x18	75004122	-	2	3	-	4	

Tabulka 40: Nabízené univerzální odkládací tácy pro základny Solaris 2000

Pro základny Solaris 4000

	Č. výrobku Univerzální		zální	Dvoupatrová
		18 x 30	36 x 24	18 x 30
Univerzální odkládací tác 11x14 / 12x14	75004123	-	-	-
Univerzální odkládací tác 9x18	75004122	3	6	6

Tabulka 41: Nabízené univerzální odkládací tácy pro základny Solaris 4000

Pro základny Solaris 2000 I / 2000 R

	Č. výrobku	Univerzální
		11 x 14
Univerzální odkládací tác 11x14 / 12x14	75004123	1
Univerzální odkládací tác 9x18	75004122	-

Tabulka 42: Nabízené univerzální odkládací tácy pro základny Solaris 2000 I / 2000 R

Pro základny Solaris 4000 I / 4000 R

	Č. výrobku	Univerzální
		18 x 18
Univerzální odkládací tác 11x14 / 12x14	75004123	-
Univerzální odkládací tác 9x18	75004122	2

Tabulka 43: Nabízené univerzální odkládací tácy pro základny Solaris 4000 I / 4000 R

Příslušenství pro univerzální odkládací tácy

	Č. výrobku	Univerzální
Náhradní O-kroužky (univerzální odkládací tác)	75004132	-

Tabulka 44: Nabízené příslušenství pro univerzální odkládací tácy

1. 2. 11. Adhezní podložky

Obecné pokyny k výběru adhezních podložek:

- Podložky o velikosti 170x280 mm jsou optimálně vhodné pro základny 11x14 a 12x14.
- Podložky o velikosti 200x200 mm jsou optimálně vhodné pro větší základny 18x18, 18x24, 30x18 a 36x24.
- Všechny podložky je možné přiříznout na jakoukoli požadovanou velikost.

Pro základny Solaris 2000

	Č. výrobku	Univerzální		í	Dvoupatrová univerzální	
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Podložka s vysokou přilnavostí, 200x200 mm, sada: 2 ks	75004126	1	4	6	2	8
Podložka s vysokou přilnavostí, 280x170 mm, sada: 2 ks	75004127	2	2	4	4	4
Podložka s nízkou přilnavostí, 200x200 mm, sada: 2 ks	75004111	1	4	6	2	8
Podložka s nízkou přilnavostí, 280x170 mm, sada: 2 ks	75004117	2	2	4	4	4

Podložky jsou prodávány v sadách po 2 kusech. Počet podložek, který je vhodný pro danou základnu, se nevztahuje na velikost sady, ale na jednotlivé podložky.

Tabulka 45: Nabízené adhezní podložky pro základny Solaris 2000

Pro základny Solaris 4000

	Č. výrobku	Univerzální		Dvoupatrová
		18 x 30	36 x 24	18 x 30
Podložka s vysokou přilnavostí, 200x200 mm, sada: 2 ks	75004126	6	12	12
Podložka s vysokou přilnavostí, 280x170 mm, sada: 2 ks	75004127	6	9	12
Podložka s nízkou přilnavostí, 200x200 mm, sada: 2 ks	75004111	6	12	12
Podložka s nízkou přilnavostí, 280x170 mm, sada: 2 ks	75004117	6	9	12

Podložky jsou prodávány v sadách po 2 kusech. Počet podložek, který je vhodný pro danou základnu, se nevztahuje na velikost sady, ale na jednotlivé podložky.

Tabulka 46: Nabízené adhezní podložky pro základny Solaris 4000



Pro základny Solaris 2000 I / 2000 R

	Č. výrobku	Univerzální
		11 x 14
Podložka s vysokou přilnavostí, 200x200 mm, sada: 2 ks	75004126	1
Podložka s vysokou přilnavostí, 280x170 mm, sada: 2 ks	75004127	2
Podložka s nízkou přilnavostí, 200x200 mm, sada: 2 ks	75004111	1
Podložka s nízkou přilnavostí, 280x170 mm, sada: 2 ks	75004117	2

Podložky jsou prodávány v sadách po 2 kusech. Počet podložek, který je vhodný pro danou základnu, se nevztahuje na velikost sady, ale na jednotlivé podložky.

Tabulka 47: Nabízené adhezní podložky pro základny Solaris 2000 I / 2000 R

Pro základny Solaris 4000 I / 4000 R

	Č. výrobku	Univerzální
		18 x 18
Podložka s vysokou přilnavostí, 200x200 mm, sada: 2 ks	75004126	4
Podložka s vysokou přilnavostí, 280x170 mm, sada: 2 ks	75004127	2
Podložka s nízkou přilnavostí, 200x200 mm, sada: 2 ks	75004111	4
Podložka s nízkou přilnavostí, 280x170 mm, sada: 2 ks	75004117	2

Podložky jsou prodávány v sadách po 2 kusech. Počet podložek, který je vhodný pro danou základnu, se nevztahuje na velikost sady, ale na jednotlivé podložky.

Tabulka 48: Nabízené adhezní podložky pro základny Solaris 4000 I / 4000 R

Příslušenství

1. 2. 12. Obecné příslušenství

1	. 2. 12. Obecné příslušenství	í	No.
		Č. výrobku	- V
	Aplikační váleček pro adhezní podložky	75004124	
	Šroubovák pro díly příslušenství	75004131	
	Rozdělovač plynu (Solaris 2000 I / 2000 R)	SK2000-8GM	
	Rozdělovač plynu (Solaris 4000 I / 4000 R)	SK4000-8GM	
	Sada náhradních dílů pro držák baněk (šrouby)	SK0010	
	Sada náhradních dílů pro základnu Solaris 2000, 2000 I/R a 4000 I/R (šrouby základny, nářadí)	SK0100	
	Sada náhradních dílů pro základnu SK4000 (šrouby základny, nářadí)	SK0101	

Tabulka 49: Nabízené obecné příslušenství

1.3. Normy a směrnice

Region	Směrnice	Normy
Evropa	2006/42/ES	EN 61010-1
	Směrnice o strojních zařízeních	IEC 61010-2-051
	2011/65/EU RoHS	EN 61326-1 třída B
	Omezení používání určitých	EN ISO 14971
	nebezpečných látek v elektrických a elektronických zařízeních.	EN ISO 9001
	Ochranné cíle:	<u>Jen pro Solaris 2000 I / 2000 R / 4000 I /</u> 4000 R: IEC 61010-2-010
	2014/35/EU o nízkonapěťových zařízeních	
	2014/30/EC Elektromagnetická kompatibilita (EMC)	
Severní Amerika		ANSI/UL 61010-1
		IEC 61010-2-051
		IEC 61326-1 třída B
		CFR 47 FCC 15 EMV
		EN ISO 14971
		EN ISO 9001
		<u>Jen pro Solaris 2000 I / 2000 R / 4000 I /</u> <u>4000 R:</u> IEC 61010-2-010
Japonsko		IEC 61010-1
		IEC 61010-2-051
		IEC 61326-1 třída B
		EN ISO 14971
		EN ISO 9001
		Jen pro Solaris 2000 I / 2000 R / 4000 I /
		4000 R: IEC 61010-2-010

Tabulka 50: Normy a směrnice

UPOZORNĚNÍ: Toto zařízení bylo testováno a splňuje mezní hodnoty pro digitální zařízení třídy A podle předpisů FCC, část 15. Tyto mezní hodnoty jsou navrženy tak, aby zajišťovaly přiměřenou ochranu proti nežádoucímu rušení při použití zařízení v komerčním prostředí. Toto zařízení generuje, používá a může vyzařovat rádiové vlnění. Není-li instalováno a používáno podle pokynů výrobce, může způsobit nežádoucí rušení rádiové komunikace. Použití tohoto zařízení v obytném prostředí může způsobit nežádoucí rušení. V takovém případě je uživatel povinen rušení na své náklady odstranit.

1.4. Přehled výrobku

1. 4. 1. Solaris 2000



① Montážní deska základny; ② Grafické uživatelské rozhraní; ③ Síťový vypínač; ④ Síťový konektor;
 ⑤ Pojistka; ⑥ Ethernetový konektor; ⑦ USB konektor

Obraz 1: Přehled Solaris 2000

1. 4. 2. Solaris 4000



① Montážní deska základny; ② Grafické uživatelské rozhraní; ③ Síťový vypínač; ④ Síťový konektor;
 ⑤ Pojistka (s resetem); ⑥ Ethernetový konektor; ⑦ USB konektor

Obraz 2: Přehled Solaris 4000

Přehled výrobku

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



① Montážní deska základny; ② Grafické uživatelské rozhraní; ③ Síťový vypínač; ④ Kabelová průchodka;
 ⑤ Síťový konektor; ⑥ Pojistka (s resetem); ⑦ Ethernetový konektor; ⑧ USB konektor; ⑨ Kryt

Obraz 3: Přehled Solaris 2000 I / 2000 R

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



① Montážní deska základny; ② Grafické uživatelské rozhraní; ③ Síťový vypínač; ④ Kabelová průchodka;
 ⑤ síťový konektor; ⑥ Pojistka (s resetem); ⑦ Ethernetový konektor; ⑧ USB konektor; ⑨ Kryt

Obraz 4: Přehled Solaris 4000 I / 4000 R

1.4.5. Přípojky

Otevřená míchadla



① Síťový konektor; ② Pojistka; ③ Ethernetový konektor

Obraz 5: Zadní pohled (vlevo: Solaris 4000, vpravo: Solaris 2000)

Inkubovaná míchadla (jen vyhřívaná) a chlazená míchadla (chlazená a vyhřívaná)



1Síťový konektor; 2 Pojistka; 3 Ethernetový konektor; 4 Odtoková hadice

Obraz 6: Zadní pohled (vlevo: Solaris 2000 I / 2000 R, vpravo: Solaris 4000 I / 4000 R)

Síťový konektor

Míchadlo vyžaduje napájecí zdroj, který odpovídá jeho specifikaci. Síťové kabely jsou součástí dodávky.

Č. výrobku	Míchadlo	Technické údaje
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

Obraz 7: Specifikace napájecího zdroje míchadel

VÝSTRAHA Poškození v důsledku připojení k nesprávnému napájecímu zdroji nebo použití nesprávného konektoru. Zkontrolujte, že je míchadlo připojeno jen do elektrické zásuvky řádně uzemněné podle platných norem pro elektrickou bezpečnost. Neprovozujte míchadlo s poškozeným nebo nedostatečně dimenzovaným síťovým kabelem.

Síťová zástrčka musí být vždy volně přístupná.

V rámci ochrany před zásahem elektrickým proudem zkontrolujte, že je plocha kolem míchadla suchá. Dojde-li nedopatřením k rozlití nebo vystříknutí kapalin, odpojte míchadlo od elektrické sítě, odstraňte rozlitou kapalinu a odstraňte všechna zdravotní nebo biologická nebezpečí, až pak pokračujte v práci.

Při delším nepoužívání odpojte míchadlo od elektrické sítě.

<u>Pojistka</u>

Pojistku je po její aktivaci nadměrným proudem možné resetovat. Míchadlo je možné uvést znovu do provozu až po resetu pojistky do správné polohy.

Ethernet

Míchadlo je vybaveno ethernetovým konektorem RJ45, který je možné použít k připojení do lokální datové sítě (LAN). Do ethernetového konektoru RJ45 je dovoleno připojovat jen zařízení, která splňují normu IEC 60950-1. Přes ethernetový konektor RJ45 je později možné nahrát dostupné softwarové aktualizace.

Kabelové průchodky

Míchadla s regulací teploty mají po stranách dva otvory, které označujeme jako kabelové průchodky. Tyto kabelové průchodky jsou zpravidla uzavřené krytkami. Po sejmutí krytky můžete otvorem zavést další zařízení, např. přídavný teplotní senzor, nebo protáhnout kabely a hadice, např. hadice rozdělovače plynu.

<u>USB</u>

Míchadlo je vybaveno dvěma konektory USB-A 2.0, které je možné použít k připojení USB flash disku. Do USB konektoru je dovoleno připojovat jen zařízení, která splňují normu IEC 60950-1.

Interní ochrana proti přehřátí

Míchadla s regulací teploty jsou vybavena termostatem s ručním resetem, který je umístěný vedle topných prvků. Toto termické zařízení vypne v případě výpadku cirkulačních ventilátorů ohřev, aby uvnitř míchadla nedošlo k přílišnému nárůstu teploty.

Aktivovaný termostat je před dalším provozem míchadla nutné ručně resetovat. Tento krok může provést jen zákaznický servis společnosti Thermo Fisher Scientific.

2. Přeprava a instalace



Za řádnou instalaci míchadla odpovídá sám zákazník.

UPOZORNĚNÍ

Ihned při dodání zkontrolujte přepravní obal. Před vybalením výrobku pečlivě zkontrolujte, zda během přepravy nedošlo k žádnému poškození. Zjištěná poškození nechte zaznamenat a podepsat dopravcem do vaší kopie dodacího listu.

Opatrně otevřete obal a zkontrolujte, že jsou k dispozici všechny komponenty ("Tabulka 51: Obsah dodávky"), až poté obalový materiál zlikvidujte. Zjistíte-li po rozbalení poškození výrobku, oznamte tuto skutečnost dopravci a požádejte o přezkoumání škody.

Důležité: Nepožádáte-li do několika dní po obdržení zásilky o přezkoumání škody, je dopravce zbaven odpovědnosti za vzniklé škody. O přezkoumání škody musíte požádat.

2.1. Vybalení

Při vybalení zkontrolujte podle seznamu balení, že jste obdrželi kompletní obsah dodávky. Obalový materiál nelikvidujte, dokud jste se nepřesvědčili o kompletnosti všech dílů.

Obsah dodávky

Artikl	Množství
Míchadlo	1
Napájecí kabel	1
Univerzální základna	1*
Šrouby základny	
- Solaris 4000	4*
- ostatní provedení	3*
Návody (výtisk, cz)	1
Návod k použití na USB flash disku	1
Upevňovací nářadí (klíč s T-rukojetí)	1
* U míchadel 2000 I, 2000 R, 4000 I a 4000 R je už základna předinstalovaná v míchadle. Základna ani šrouby nejsou k dispozici jako samostatné díly.	

Tabulka 51: Obsah dodávky

Nebyly-li dodány všechny díly, kontaktujte společnost Thermo Fisher Scientific.

2. 2. Místo instalace

POZOR V důsledku degradace plastů vystavených působení ultrafialového záření může dojít ke snížení bezpečnosti zařízení. Chraňte míchadlo a díly příslušenství z plastu před přímým slunečním zářením a ostatními zdroji ultrafialového záření.

Postavte míchadlo na rovný stůl nebo lavici, které unesou hmotnost běžícího míchadla s příslušenstvím a vzorky. Umístěte míchadlo v blízkosti elektrické zásuvky, která splňuje požadavky typového štítku. Ponechte okolo zařízení dostatečný prostor pro volné proudění vzduchu, instalované části příslušenství a vyšší komfort obsluhy.

Při instalaci zařízení dodržte následující požadavky:

- Pro zajištění dostatečné vzdálenosti od pohyblivých dílů a pro zabránění přiskřípnutí/ pohmoždění částí těla nebo poškození okolních zařízení je u otevřených míchadel nutné ponechat okolo základny odstup 8 cm (3 palce) do všech čtyř stran. Základna otevřeného míchadla může přesahovat přes plášť míchadla.
- Míchadla způsobují vibrace. V bezpečnostní zóně se nesmí nacházet žádná citlivá zařízení, nesmí se v ní skladovat nebezpečné předměty a látky.



Obraz 8: Instalace otevřených míchadel s bezpečnostním odstupem 8 cm (3 palce)

Pro zajištění řádného odvětrání a optimálního výkonu je u míchadel s regulací teploty nutné dodržet následující odstupy:

- Inkubovaná míchadla (jen vyhřívaná) Solaris 2000 I a 4000 I vyžadují okolo pláště bezpečnostní odstup 8 cm (3 palce) do všech čtyř stran, viz Obraz 9.
- Chlazená míchadla (chlazená a vyhřívaná) Solaris 2000 R a 4000 R je možné provozovat s odstupem 8 cm (3 palce) do všech čtyř stran okolo pláště, je-li nad krytem volný prostor, jak je znázorněno na levé straně Obraz 9.

Je-li volný prostor nad chlazeným míchadlem (chlazeným a vyhřívaným) Solaris 2000 R a 4000 R zastavěný, je odstup okolo pláště nutné zvětšit do všech čtyř stran na 30 cm (12 palců), jak je znázorněno na pravé straně Obraz 9.



Obraz 9: Instalační vzdálenosti pro inkubovaná míchadla (jen vyhřívaná) a chlazená míchadla (chlazená a vyhřívaná) s volným prostorem nad zařízením (vlevo), a pro zastavěná chlazená míchadla (chlazená a vyhřívaná) (vpravo)

VÝSTRAHA Během míchání se nesmí v této oblasti nacházet žádné nebezpečné látky.

Podklad musí:

- » být čistý,
- » být stabilní, robustní, tvrdý a bez vibrací,
- » umožňovat vodorovné ustavení míchadla,
- » unést hmotnost míchadla.
- Míchadlo je dovoleno používat jen ve vnitřních prostorech.
- Míchadlo nesmí být vystaveno teplu ani silnému slunečnímu záření.
- Místo instalace musí být neustále dobře odvětráváno.
- Síťová zástrčka musí být vždy volně přístupná.

2.3. Přeprava

POZOR Zranění v důsledku pádu míchadla. Míchadlo zvedejte vždy z obou stran. Míchadlo nezvedejte za čelní kryt ani za instalovanou základnu. Míchadlo s regulací teploty nikdy nezvedejte za držadlo průhledného krytu.

POZOR Bezpečný provoz míchadla ohrožuje kondenzace vody v plášti. K ní dochází při přemístění míchadla z chladnějšího místa na teplejší místo s vlhčím vzduchem, např. ze skladu nebo přepravního vozu do laboratoře. Nechte zařízení před uvedením do provozu alespoň 2 hodiny vyschnout.

UPOZORNĚNÍ Před manipulací s otevřenými míchadly vždy demontujte základnu, sejměte náklad a všechny díly příslušenství. Neodstraníte-li základnu, může případně dojít k poškození montážní desky základny nebo míchacího mechanismu. U míchadel s regulací teploty stačí ze základny sejmout náklad.

UPOZORNĚNÍ Při nárazu se může zařízení poškodit. Míchadlo přepravujte ve vzpřímené poloze a, je-li to možné, v originálním obalu.

Manipulace s míchadlem

Při manipulaci s míchadlem vždy dodržte tyto pokyny:

- Míchadlo zvedejte vždy z obou stran, nikdy však zepředu a zezadu.
- Míchadlo musí s ohledem na jeho hmotnost přepravovat dostatečný počet osob, vždy však alespoň dvě osoby (další informace najdete v oddíle "Technické specifikace" na straně 11).







Obraz 10: Přenášení míchadla

VÝSTRAHA Míchadlo zvedejte vždy z obou stran. Míchadlo nezvedejte za čelní ani zadní stranu. Míchadla jsou těžká (další informace najdete v oddíle "Technické specifikace" na straně 11). Zvednutí a přenos míchadla vyžaduje spolupráci alespoň 2 osob.

POZOR Míchadlo s regulací teploty nikdy nezvedejte a nepřenášejte za držadlo průhledného krytu. Držadlo se může ulomit a míchadlo může spadnout na zem a způsobit těžká zranění osob.

Nivelace

2.4. Nivelace

Míchadlo samotné není vybaveno žádným nivelačním přípravkem. Řádná instalace míchadla vyžaduje vodorovný podklad.

UPOZORNĚNÍ Mezi podklad a nohy míchadla nevkládejte žádné distanční podložky ani jiné ploché předměty.

2.5. Síťový konektor

Míchadlo vyžaduje napájecí zdroj, který odpovídá jeho specifikaci. Síťové kabely jsou součástí dodávky. Další informace najdete na Obraz 7 na straně 47.

VÝSTRAHA Poškození v důsledku připojení k nesprávnému napájecímu zdroji nebo použití nesprávného konektoru. Zkontrolujte, že je míchadlo připojeno jen do řádně uzemněné elektrické zásuvky. Neprovozujte míchadlo s poškozeným nebo nedostatečně dimenzovaným síťovým kabelem.

UPOZORNĚNÍ Elektromagnetické záření může rušit zobrazení na displeji. Nedochází však k poškození ani omezení/změně funkce zařízení. Aby nedocházelo k rušení elektromagnetickým zářením, nepoužívejte v bezprostřední blízkosti zařízení žádné mobilní přístroje, např. mobilní telefony. Neprovozujte zařízení v jednom proudovém okruhu s jinými zařízeními s vysokým elektrickým výkonem. Neprovozujte více zařízení připojených do společné zásuvkové lišty.

Míchadlo připojte k elektrické síti následujícím způsobem:

- 1. Vypněte síťový vypínač na pravé straně.
- 2. Zkontrolujte, že kabel splňuje bezpečnostní předpisy vaší země.
- 3. Zkontrolujte, že síťové napětí a frekvence souhlasí s údaji na typovém štítku.

Síťová zástrčka musí být vždy volně přístupná.

V rámci ochrany před zásahem elektrickým proudem zkontrolujte, že je plocha kolem míchadla suchá. Dojde-li nedopatřením k rozlití nebo vystříknutí kapalin, odpojte míchadlo od elektrické sítě a odstraňte rozlité kapaliny, až pak pokračujte v práci.

Při delším nepoužívání odpojte míchadlo od elektrické sítě.

2.6. Základní nastavení

Před použitím zařízení proveďte následující kroky:

1. Zapněte zařízení. Zobrazí se logo Thermo Scientific. Stiskněte tlačítko Start Setup.



Obraz 11: Výzva k provedení základního nastavení

2. Zvolte požadovaný jazyk na obrazovce Jazyk. Stiskněte tlačítko Pokračovat.



Obraz 12: Základní nastavení - Jazyk

 V dialogovém okně Název zařízení můžete případně zadat název zařízení. Stiskněte tlačítko Pokračovat.

	Unit Nar	ne	
Unit Name:	Tap to enter		
E	Back	Next	

Obraz 13: Základní nastavení - Název zařízení

 Do textového pole můžete případně zadat město a stát. Po zadání prvních třech znaků se v textovém poli zobrazí seznam vhodných návrhů. Zvolte případně některý z návrhů a stiskněte tlačítko Pokračovat.

Set Re	gion
Enter the city and cou	intry closest to your time zone
Tap to enter	
	Set Re

Obraz 14: Základní nastavení - Region

5. Zvolte požadovaný formát data. Stiskněte tlačítko Pokračovat.

Set Date Dis	splay Format
MMI	M/DD/YYYY
	MMM/YYYY
O YYY	Y/MMM/DD
Back	Next

Obraz 15: Základní nastavení - Nastavení formátu data

 Pomocí jednotlivých segmentů nastavovacího kolečka nastavte aktuální datum. Stiskněte tlačítko Pokračovat.



Obraz 16: Základní nastavení - Nastavení data

 Zvolte požadovaný formát času a pomocí jednotlivých segmentů nastavovacího kolečka nastavte aktuální čas. Stiskněte tlačítko Pokračovat.



Obraz 17: Základní nastavení - Nastavení času

 Jen inkubovaná míchadla (jen vyhřívaná) a chlazená míchadla (chlazená a vyhřívaná): zvolte požadovanou jednotku teploty. Stiskněte tlačítko Pokračovat.

Temper	ature Units
	°C
C) *F
Back	Next

Obraz 18: Základní nastavení - Volba jednotky teploty

 Jen inkubovaná míchadla (jen vyhřívaná) a chlazená míchadla (chlazená a vyhřívaná): Stiskněte modrou šipku a nastavte alarm nadměrné teploty a mezní hodnoty pro nouzové vypnutí. Stiskněte tlačítko Pokračovat.



Obraz 19: Základní nastavení - Nastavení mezních hodnot alarmu příliš vysoké a příliš nízké teploty

10. Zopakujte postup pro mezní hodnoty alarmu příliš nízké teploty. Stiskněte tlačítko Pokračovat.

 Chcete-li pokračovat bez nastavení přístupového hesla administrátora, nechte nastavenou volbu Bez řízení přístupu a stiskněte tlačítko Pokračovat.

Chcete-li nastavit přístupové heslo administrátora, stiskněte volbu S řízením přístupu.

	Acces	s Mode
•	Open Mode Select to allow full access to	c all functionality for all users.
0	Secure Mode Select to require users to er changes to stored paramete (All users still able to use pro	nter the admin passcode to make ers, programs, and settings. grams and set new run parameters.
1	Pack	

Obraz 20: Základní nastavení - Volba režimu přístupu

 V zobrazené výzvě k zadání přístupového hesla zadejte pomocí klávesnice přístupové heslo administrátora a stiskněte tlačítko Pokračovat.

Admin	Passcode
The unit has shipped with a s	tandard passcode. If you would
like to change the standard adr	min passcode, please enter a new
passco	ide below.
If you do not wish to change th	e passcode at this time, leave the
fields below blank ar	nd tap the Next button.
Back	Next

Obraz 21: Základní nastavení - Volba režimu přístupu

- 13. V další výzvě k zadání přístupového hesla zadejte své přístupové heslo administrátora.
- Ve třetí výzvě k zadání přístupového hesla zadejte pro potvrzení ještě jednou své přístupové heslo administrátora.

Admin Passcode			Admin Passcode							
If you do not wis fields	sh to change below blank	the pa and ta	sscode a ap the Ne	at this time, leave the ext button.	F	Please re-ent	er the n	ew passcode		
*	*	*		* *	*	*	*	*	*	
	7	8	9			7	8	9		
	4	5	6			4	5	6		
	1	2	3			1	2	3		
	0	Cl	ear			0	Cle	ear		

Obraz 22: Základní nastavení - Zadání a potvrzení nového přístupového hesla administrátora

Základní nastavení

15. Na zobrazené potvrzovací obrazovce stiskněte tlačítko Pokračovat.

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

Obraz 23: Základní nastavení - Potvrzení změny přístupového hesla administrátora

 Zobrazí se obrazovka s obecnými fyzickými instalačními pokyny k míchadlu. Stiskněte tlačítko Pokračovat.

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					

Obraz 24: Základní nastavení - Instalace a připojení

 Zobrazí se obrazovka s informací o dokončené konfiguraci. Stiskněte tlačítko Pokračovat, proces se dokončí.

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

Obraz 25: Základní nastavení - Dokončení kalibrace nebo základního nastavení

UPOZORNĚNÍ U inkubovaných míchadel (jen vyhřívaných) a chlazených míchadel (chlazených a vyhřívaných) můžete stisknout tlačítko **Kalibrovat** a provést proces popsaný v oddíle "4. 6. Kalibrace teploty" na straně 126.

2.7. Skladování



POZOR

Před uskladněním míchadla a příslušenství celý systém vyčistěte a v případě nutnosti proveďte jeho dezinfekci nebo dekontaminaci. Míchadlo a příslušenství neuskladňujte, dokud není jednoznačně definovaný stav kontaminace. Nejste-li si jisti postupem, kontaktujte zákaznický servis společnosti Thermo Fisher Scientific ("Čištění" na straně 123, "Dezinfekce" na straně 124 a "Dekontaminace" na straně 124).

- Před uskladněním míchadla a příslušenství vše vyčistěte a v případě nutnosti proveďte dezinfekci a dekontaminaci.
- Před uskladněním musí být míchadlo i příslušenství zcela suché.
- Uskladněte míchadlo na čistém, bezprašném místě.
- Uskladněte míchadlo tak, aby stálo na nohách.
- Uskladněte míchadlo na místě bez přímého slunečního záření.

2.8. Odeslání



Před odesláním míchadla a příslušenství celý systém vyčistěte a v případě nutnosti proveďte jeho dezinfekci nebo dekontaminaci. Míchadlo a příslušenství neuskladňujte, dokud není jednoznačně definovaný stav kontaminace. Nejste-li si jisti postupem, kontaktujte zákaznický servis společnosti Thermo Fisher Scientific ("Čištění" na straně 123, "Dezinfekce" na straně 124 a "Dekontaminace" na straně 124).

Před odesláním míchadla dodržte následující:

- Míchadlo musí být vyčištěné a dekontaminované.
- Provedení dekontaminace musí být potvrzeno dekontaminačním certifikátem.
 Dekontaminační certifikát si můžete vyžádat u zákaznického servisu společnosti Thermo Fisher Scientific.

3. Obsluha

3.1. Zapnutí/vypnutí

Stiskněte síťový vypínač na pravé straně, míchadlo se zapne, resp. vypne ((I) nebo (0)).

Během bootování se zobrazuje logo Thermo Scientific.

V připraveném stavu se na ovládacím panelu zobrazuje aktuální provozní stav míchadla.

3. 2. Grafické uživatelské rozhraní

Hlavní obrazovka je standardní obrazovkou grafického uživatelského rozhraní míchadla. Představuje výchozí bod pro provoz míchadla.

Na hlavní obrazovce můžete:

- nastavit nejdůležitější provozní parametry, např. otáčky míchadla, délku míchání a teplotu, (jen míchadla s regulací teploty)
- spustit a zastavit míchadlo,
- zobrazit informace k provoznímu stavu a zpracovat alarmová a výstražná hlášení,
- přejít na další obrazovky s informacemi k provoznímu stavu a s možnostmi nastavení.

Obsah obrazovky otevřených míchadel a míchadel s regulací teploty se mírně liší.

Na Obraz 26 vidíte příklad hlavní obrazovky otevřeného míchadla.



Obraz 26: Grafické uživatelské rozhraní - Hlavní obrazovka otevřeného míchadla

Hlavní obrazovka míchadla s regulací teploty obsahuje také prvky pro regulaci teploty, jak je znázorněno na Obraz 27.



Obraz 27: Grafické uživatelské rozhraní - Hlavní obrazovka míchadla s regulací teploty

3. 2. 1. Nastavení nejdůležitějších provozních parametrů míchadla

Z hlavní obrazovky přejdete jedním stisknutím k nastavovacím obrazovkám nejdůležitějších provozních parametrů míchadla.

Nastavení otáček

 Stiskněte pole Skutečné otáčky (© na Obraz 26 nebo ⑦ na Obraz 27), otevře se obrazovka Požadované otáčky, která je znázorněná na Obraz 28.



Obraz 28: Požadované otáčky

 Stiskněte šipky nad nebo pod každým segmentem nastavovacího kolečka a nastavte požadované otáčky. Alternativně je požadované otáčky možné nastavit i otáčením jednotlivých segmentů nastavovacího kolečka.

3. Stiskněte tlačítko Uložit.

UPOZORNĚNÍ Zobrazí-li se vyskakovací okno s informací, že požadované otáčky leží mimo přípustný rozsah, je vaše nastavení mimo rozsah provozních otáček, které míchadlo podporuje, viz oddíl "Technické údaje" na straně 11. Opravte nastavení a pokračujte.

4. Potvrďte následující dialogové okno Nastavení byla uložena, vrátíte se na hlavní obrazovku.

Výpočet kruhové dráhy: Výpočet přibližných otáček z nastavení předchozího zařízení

V dialogovém okně Požadované otáčky můžete také stisknout tlačítko **Výpočet otáček**, otevře se obrazovka **Výpočet kruhové dráhy**. Tuto funkci je možné využít při přechodu z míchadla s jinou mechanickou nástavbou na některé míchadlo konstrukční řady Solaris popsané v tomto návodu. Tato funkce zohledňuje rozdíl velikostí kruhové dráhy a nabídne hrubý odhad požadovaných otáček, které jsou nutné pro dosažení podobného výsledku.

UPOZORNĚNÍ Výsledný přepočet ihned nepoužívejte pro výrobní vzorky, vždy proveďte nejprve několik ověřovacích testů. V závislosti na aktuálním smykovém namáhání mohou buňky růst rychleji nebo pomaleji, vytvářet jiné proteiny, poškodit se (odumřít), protože při přechodu na jinou kruhovou dráhu se mění fyzikální podmínky.

1. Zvolte měrné jednotky (milimetry nebo palce).



Obraz 29: Výpočet kruhové dráhy

 Zvolte velikost předchozí kruhové dráhy. Stiskněte tlačítko Vypočítat. Pro výpočet otáček se zobrazí následující okno.



Obraz 30: Vypočtená kruhová dráha

- 3. Stiskněte tlačítko OK.
- Doporučené otáčky je možné vypočítat pro různě velikou kruhovou dráhu dle potřeby zákazníka. Pro výpočet vlastní kruhové dráhy zvolte možnost Jiná.
- 5. Zadejte velikost kruhové dráhy.



Obraz 31: Výpočet vlastní kruhové dráhy

6. Stiskněte tlačítko Vypočítat.

Zobrazí se okno Výpočet otáček a ukáže otáčky pro zadanou velikost kruhové dráhy.

Nastavení procesního času

Míchadlo je možné provozovat v trvalém provozu nebo dle nastaveného času. V trvalém provozu je míchadlo možné zastavit podle potřeby ručně. V provozu dle nastaveného času se pohon míchadla zastaví automaticky po uplynutí času nastaveného na časovači. Časovač je možné zobrazit následujícím způsobem:

- uplynulý čas: jak dlouho už míchadlo běží od okamžiku, kdy jste stiskli tlačítko Start, nebo
- zbývající čas: jak dlouho ještě míchadlo poběží, než uplyne čas nastavený na časovači.
- Stiskněte pole Čas (⑤ na Obraz 26 a Obraz 27), otevře se obrazovka Časový režim, která je znázorněná na Obraz 32.
- 2. Zvolte provoz dle nastaveného času.



Obraz 32: Volba časového režimu

- 3. Zadejte v Nastavení času hodiny a minuty.
- 4. Stiskněte tlačítko Uložit.
- 5. Pro návrat na hlavní obrazovku potvrďte následující dialogové okno Nastavení byla uložena.

Nastavení teploty

U míchadel s regulací teploty můžete v poli Požadovaná teplota kdykoli upravit teplotu aplikace.

- Stiskněte pole Požadovaná teplota (spodní oblast poz. 6 na Obraz 27), otevře se obrazovka Požadovaná teplota, která je znázorněná na Obraz 33.
- Stiskněte šipky nad nebo pod každým segmentem nastavovacího kolečka a nastavte standardní teplotu, kterou má míchadlo udržovat. Alternativně je požadovanou teplotu možné nastavit i otáčením jednotlivých segmentů nastavovacího kolečka.



Obraz 33: Nastavení požadované teploty: obrazovka s kalibračním offsetem

3. Stiskněte tlačítko Uložit.

UPOZORNĚNÍ Zobrazí-li se vyskakovací okno s informací, že požadovaná teplota leží mimo přípustný rozsah, je vaše nastavení mimo rozsah provozních teplot, které míchadlo podporuje, viz oddíl "Technické údaje" na straně 11. Opravte nastavení a pokračujte.

UPOZORNĚNÍ Zobrazí-li se vyskakovací okno s informací, že zvolená teplota leží vzhledem k aktuální teplotě okolí mimo přípustný rozsah, je teplota okolí mimo rozsah provozních teplot míchadla, viz oddíl "Technické údaje" na straně 11. Stiskněte tlačítko OK pro potvrzení, že jste informaci vzali na vědomí a chcete pokračovat, nebo zvolte jinou teplotu.

UPOZORNĚNÍ Obrazovka Požadovaná teplota zobrazuje kalibrovanou teplotu a hodnotu offsetu nastavenou během kalibrace, jak je znázorněno v levé spodní části na Obraz 33. Kalibrační proces je popsaný v oddíle "Kalibrace teploty" na straně 126.

4. Pro návrat na hlavní obrazovku potvrďte následující dialogové okno Nastavení byla uložena.

UPOZORNĚNÍ U míchadel s regulací teploty dochází před dosažením požadované teploty k tzv. překmitu. Překmit teploty znamená, že teplota komory nejprve dosáhne o trochu vyšší (nebo nižší) teploty a pak se přibližuje požadované teplotě. V této přibližovací fázi se příslušný překmit na displeji nezobrazuje. Místo toho se na displeji zobrazuje až do dosažení požadované teploty průběžný nárůst (nebo pokles) teploty komory.

Spuštění a zastavení míchadla

1. Pro spuštění míchadla stiskněte tlačítko Start.

Tlačítko Start se změní na tlačítko Pauza.



Obraz 34: Tlačítka Start resp. Stop

2. Pro zastavení míchadla stiskněte tlačítko Stop.

3. 2. 2. Provozní stav

Nachází-li se míchadlo v řádném provozním stavu, zobrazí se na ovládacím panelu v oblasti



Provozní stav a statistické údaje (⑦ na Obraz 26 a ⑧ na Obraz 27) zelený symbol srdce. Stisknutím zeleného symbolu srdce se otevře obrazovka Provozní stav. Obrazovka Provozní stav obsahuje statistické údaje k provozu míchadla: míchací

a provozní hodiny celkem a míchací a provozní hodiny poslední relace. U míchadel s regulací teploty se zobrazují i hodiny chladicího a topného procesu.

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

Obraz 35: Statistické údaje míchadla: otevřené míchadlo (vlevo) a míchadlo s regulací teploty (vpravo)

Výstražné hlášení

Při vydání výstražného hlášení se na ovládacím panelu v horní části aktuální obrazovky zobrazí žlutý pruh s běžícím textem. Zároveň zazní akustický signál. Žlutý výstražný pruh

s běžícím textem zmizí po druhém průběhu hlášení. Skutečnost, že je aktivní jedno nebo více alarmových hlášení míchadla, indikuje jen žlutý výstražný trojúhelník.



Součástí výstražného trojúhelníku je modrý kroužek s bílým okrajem, ve kterém se zobrazuje počet aktivních výstražných hlášení. Stisknutím výstražného trojúhelníku v oblasti Info a provozní stav (⑦ na Obraz 26 a ⑧ na Obraz 27) se otevře obrazovka se seznamem všech aktuálně aktivních výstražných hlášení. Nejnovější výstražné hlášení se zobrazí v rozšířeném režimu, takže vidíte všechny informace, jak je znázorněno na Obraz 36. Seznamem můžete rolovat, klepnutím na libovolný záznam otevřete rozšířené zobrazení se všemi informacemi.



Obraz 36: Seznam výstražných hlášení

Aktivní výstražné hlášení můžete vybrat stisknutím zaškrtávacího pole vedle položky výstražného hlášení. Po stisknutí tlačítka **Potvrdit** se systém pokusí výstražné hlášení ze seznamu smazat. Po smazání všech výstražných hlášení se symbol provozního stavu opět změní na původní symbol zeleného srdce.

Alarmové hlášení

Při vydání alarmového hlášení se míchadlo okamžitě zastaví, aby nedošlo k poškození vzorků



a/nebo samotného zařízení. Před opětovným spuštěním je nutné potvrdit alarmové hlášení na ovládacím panelu.

Při vydání alarmového hlášení se na ovládacím panelu v horní části aktuální obrazovky zobrazí červený pruh. V oblasti Info a provozní stav (⑦ na Obraz 26 a ⑧ na Obraz 27) se zobrazí červený zvonek se zvukovými vlnami do obou stran. Zároveň zazní konstantní akustický signál.

Pod červeným pruhem běží textová zpráva se souhrnnou informací k aktuálnímu alarmovému hlášení. Na pravé straně se zobrazí tlačítko Ztlumení, jehož stisknutím je možné akustický signál dočasně vypnout. Není-li příčina alarmu během stanoveného času ztlumení alarmu odstraněna, zazní akustický signál znovu. Čas ztlumení alarmu je možné stanovit v nastaveních. Další informace najdete v oddíle "Alarmová/výstražná hlášení" na straně 69.

Po stisknutí zvonku v oblasti Info a provozní stav (⑦ na Obraz 26 a ⑧ na Obraz 27) se zobrazí obrazovka se všemi podrobnostmi aktuálního alarmového hlášení, jak je znázorněno na Obraz 37. Seznamem můžete rolovat, klepnutím na libovolný záznam otevřete rozšířené zobrazení se všemi informacemi.



Obraz 37: Seznam alarmových hlášení

Aktivní alarmové hlášení můžete vybrat stisknutím zaškrtávacího pole vedle položky alarmového hlášení. Stisknutím tlačítka **Potvrdit** po odstranění příčiny se systém pokusí alarmové hlášení ze seznamu smazat. Po smazání všech alarmových hlášení se symbol provozního stavu opět změní na původní symbol zeleného srdce.

<u>Chyba</u>

Dojde-li k chybě, zobrazí míchadlo chybové hlášení a okamžitě se zastaví, aby nedošlo k poškození vzorků a/nebo samotného zařízení. Obrazovka zčervená a není možná žádná další interakce. Zobrazí se chybové hlášení s chybovým kódem, jak je vidět na příkladu Obraz 38.



Obraz 38: Obrazovka s chybovým hlášením

Odstraňte problém podle následujícího popisu a pokuste se zařízení uvést opět do provozu:

- 1. Poznamenejte si chybový kód uvedený na obrazovce s chybovým hlášením.
- 2. Stiskněte tlačítko Potvrdit, akustický signál se vypne.
- 3. Restartujte míchadlo jeho vypnutím a opětovným zapnutím.
- Zobrazí-li se chybové hlášení znovu, kontaktujte zákaznický servis společnosti Thermo Fisher Scientific a sdělte chybový kód uvedený na obrazovce s chybovým hlášením.

3. 2. 3. Nastavení

Druhý symbol na navigační liště je symbol Nastavení. Po stisknutí symbolu **Nastavení** se zobrazí níže uvedená obrazovka. Obrazovka Nastavení obsahuje více tlačítek, než jich je na displeji možné najednou zobrazit. Pro zobrazení zbývajících tlačítek je proto nutné rolovat obrazovkou, jak je znázorněno na pravé straně Obraz 39.





Obraz 39: Obrazovka Nastavení

Alarmová/výstražná hlášení

Na obrazovce Alarmová/výstražná hlášení můžete určit, jak a kdy se budou vydávat alarmová a výstražná hlášení.



Obraz 40: Nastavení preferencí alarmových a výstražných hlášení

Hlasitost: Tímto jezdcem můžete upravit hlasitost akustického signálu, který zazní s alarmovým nebo výstražným hlášením. Posuňte jezdec doleva pro snížení hlasitosti, nebo doprava pro zvýšení hlasitosti, pak stiskněte tlačítko **Uložit**. Toto nastavení hlasitosti se týká jak alarmových, tak výstražných hlášení.



Obraz 41: Nastavení hlasitosti alarmových a výstražných hlášení

Alarm nadměrné teploty / alarm příliš nízké teploty (jen míchadla s regulací teploty): Pomocí těchto voleb můžete nastavit mezní hodnotu alarmu nadměrné teploty a alarmu příliš nízké teploty vzhledem k definované požadované teplotě (viz "Nastavení teploty" na straně 64). Jakmile teplota v komoře míchadla klesne pod mezní hodnotu příliš nízké teploty nebo stoupne nad mezní hodnotu nadměrné teploty, vydá míchadlo alarmové hlášení. Zvolte mezní hodnotu teploty a stiskněte tlačítko Uložit.



Obraz 42: Nastavení mezních hodnot alarmu příliš vysoké a příliš nízké teploty

UPOZORNĚNÍ Při změně požadované teploty se příslušným způsobem posunou i mezní hodnoty alarmu.

Vypnutí při nadměrné teplotě / vypnutí při příliš nízké teplotě (jen míchadla s regulací teploty): Pomocí těchto voleb můžete nastavit mezní hodnotu pro bezpečnostní vypnutí při nadměrné teplotě a při příliš nízké teplotě vzhledem k definované požadované teplotě (viz "Nastavení teploty" na straně 64). Jakmile teplota v komoře míchadla klesne pod mezní hodnotu příliš nízké teploty nebo stoupne nad mezní hodnotu nadměrné teploty, proběhne automatické vypnutí míchadla, aby nedošlo k poškození vzorků zpracovávaných v komoře. Zvolte mezní hodnotu a stiskněte tlačítko **Uložit**.

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

Obraz 43: Nastavení mezních hodnot pro bezpečnostní vypnutí při příliš vysoké a příliš nízké teplotě

UPOZORNĚNÍ Při změně požadované teploty se příslušným způsobem posunou i chybové mezní hodnoty.

Doba ztlumení: Doba ztlumení určuje, na jak dlouho se vypne akustický signál alarmového hlášení po stisknutí tlačítka Ztlumení v červeném pruhu v horní části hlavního okna (viz "Alarmové hlášení" na straně 67). Dobu ztlumení můžete nastavit na 5 minut, 10 minut a 15 minut. Standardní nastavení je 10 minut. Stiskněte tlačítko **Uložit**, provedené změny se uloží.

Snooze Timeout	5min ㅅ
	10min
	15min

Obraz 44: Nastavení doby ztlumení pro alarmová hlášení

Deaktivace kalibračních hlášení (jen míchadla s regulací teploty): Tyto volby umožňují deaktivovat oznámení, která v pravidelných intervalech připomínají nutnost provedení kalibrace měření teploty v míchadle (další informace najdete v oddíle "Kalibrace teploty" na straně 126). Můžete zvolit, zda chcete připomenutí kalibrace deaktivovat úplně, nebo jen při běžícím programu.



<u>Řízení přístupu</u>

Na obrazovce Řízení přístupu můžete přidělovat přístupová hesla administrátora, která opravňují k provádění procesů s omezenými přístupovými právy.

Míchadlo je standardně nastaveno na provoz bez řízení přístupu, tzn. že pro přístup k zařízení a jeho obsluhu není nutné zadávat přístupové heslo. Přepnutí do provozu "S řízením přístupu" vyžaduje od každého uživatele, který by chtěl měnit nastavení míchadla, přístupový kód administrátora. Je-li aktivní volba S řízením přístupu, zobrazí se v horním pravém rohu dotykové obrazovky tlačítko Přihlásit.

Následující srovnání provozu s řízením přístupu a bez řízení přístupu znázorňuje, kdy je vyžadováno zadání přístupového hesla.

Opatření	Přístupové heslo v režimu bez řízení přístupu vyžadováno	Přístupové heslo v režimu s řízením přístupu vyžadováno
Nastavení pož. otáček, doby míchání a teploty	Ne	Ne
Rozběh a zastavení míchadla	Ne	Ne
Výpočet kruhových drah	Ne	Ne
Potvrzení alarmových a výstražných hlášení	Ne	Ne
Zobrazení obrazovky Soubory a informace	Ne	Ne
Zobrazení provozního stavu	Ne	Ne
Zobrazení a export protokolu událostí a grafů	Ne	Ne
Dočasné ztlumení signálu alarmových hlášení	Ne	Ne
Použití programů	Ne	Ne
Vytváření, úprava a mazání programů	Ne	ano
Změna nastavení displeje	Ne	ano
Změna nastavení konfigurace	Ne	ano
Změna nastavení alarmů a výstrah	Ne	ano
Změna nastavení řízení přístupu	Ne	ano
Import a export programů	Ne	ano
Reset do továrního nastavení	Ne	ano
Provedení kalibrace	Ne	ano
Instalace aktualizací firmwaru	ano	ano

Tabulka 52: Závazné zadání přístupového hesla při provozu s řízením přístupu/bez řízení přístupu

Grafické uživatelské rozhraní

Obraz 46 znázorňuje obrazovku s řízením přístupu a bez řízení přístupu.

< *	Shake Lab 1 11:22:24AM Access Control		< *	Shake Lab 1 11:22:24AM Access Control	Logout Admin
¢, ∢ ∭	 Open Secure Will require passcod change certain para 	e access to meters	¢.	Open Open Secure Will require passco change certain par 	de access to ameters
	Change Admin Passcode	>		Change Admin Passcode	>
	Save	0		Save	0

Obraz 46: Řízení přístupu: Bez řízení přístupu (vlevo) a S řízením přístupu (vpravo)

Po stisknutí volby **S řízením přístupu** a následně tlačítka **Uložit** se zobrazí výzva k zadání přístupového hesla administrátora, až pak je změnu možné potvrdit. Všechna zařízení jsou dodána se stejným, továrně naprogramovaným přístupovým heslem pro administrátora. Přístupové heslo je vytištěné v návodu.

Stejným způsobem je k zadání přístupového hesla administrátora vyzván každý uživatel, který se pokusí změnit některé nastavení chráněné přístupovým heslem.

UPOZORNĚNÍ Tovární nastavení přístupového hesla administrátora je 00000.

Tlačítkem **Změnit přístupové heslo administrátora** můžete změnit přednastavené přístupové heslo. Po stisknutí tlačítka **Změnit přístupové heslo administrátora** se zobrazí výzva nejprve k zadání aktuálního přístupového hesla, pak k zadání nového přístupového hesla a nakonec k ještě jednomu zadání nového přístupového hesla pro potvrzení. Zobrazí se zpráva **Nové přístupové heslo bylo uloženo**, která informuje o úspěšném dokončení procesu.

Kalibrace

Z obrazovky Kalibrace (jen míchadla s regulací teploty) se otevírá řada obrazovek, s jejichž pomocí můžete provést kalibraci pro přesnou regulaci teploty míchadla. Tento proces je popsaný v oddíle "Kalibrace teploty" na straně 126.
Provozní parametry

Obrazovka Provozní parametry nabízí volby pro nastavení standardních provozních parametrů, se kterými se má zařízení spouštět:

Požadované otáčky: Požadované otáčky můžete nastavit na hodnotu mezi 15 a 525 ot./min. Nastavte otáčky pomocí jednotlivých segmentů nastavovacího kolečka a stiskněte tlačítko **Uložit**. Podrobné pokyny k použití obrazovky Požadované otáčky najdete v oddíle "Nastavení otáček" na straně 61.

Časový režim: Míchadlo je možné provozovat v trvalém provozu nebo dle nastaveného času. V trvalém provozu je míchadlo možné zastavit podle potřeby ručně. V provozu dle nastaveného času se pohon míchadla zastaví automaticky po uplynutí času nastaveného na časovači. Pro provoz dle nastaveného času můžete zadat standardní procesní čas a zvolit, zda se má na ovládacím panelu při běhu míchadla zobrazovat uplynulý nebo zbývající čas. Podrobné pokyny k použití obrazovky Časový režim najdete v oddíle "Nastavení procesního času" na straně 63.

Požadovaná teplota (jen míchadla s regulací teploty): Pomocí této volby můžete nastavit standardní teplotu, kterou má míchadlo při spuštění zobrazovat. Podrobné pokyny k použití obrazovky Požadovaná teplota najdete v oddíle "Nastavení teploty" na straně 64.

Automatický restart: Tato funkce znovu spustí zařízení po výpadku proudu, ke kterému došlo během normálního cyklu Start/Stop, provádění programu nebo cyklu kalibrace teploty. Je-li volba Automatický restart nastavena na **Ne**, zařízení se po výpadku proudu znovu nespustí.

<u>Displej</u>

V nastavení displeje je možné upravit různé možnosti displeje.

Jas: Pro nastavení jasu displeje použijte jezdec nebo tlačítka + / -.



Obraz 47: Nastavení jasu obrazovky

Jazyk: Pro změnu jazyka displeje stiskněte tlačítko Jazyk. Zvolte požadovaný jazyk otáčením nastavovacího kolečka a stiskněte tlačítko Uložit.



Obraz 48: Výběr jazyka obrazovky

Měrné jednotky (jen míchadla s regulací teploty): Stiskněte tlačítko Jednotky, chcete-li přepnout všechny teplotní údaje na obrazovkách grafického uživatelského rozhraní na stupně Celsia °C nebo stupně Fahrenheita °F.

Units of Measure	°C 🔨
	°F

Obraz 49: Výběr měrné jednotky teploty

Datum: Pro nastavení data stiskněte tlačítko Datum. Nastavte měsíc, den a rok pomocí jednotlivých segmentů nastavovacího kolečka a stiskněte tlačítko **Uložit**.



Obraz 50: Nastavení měsíce, data a času

Formát data: Pro změnu formátu data stiskněte tlačítko Formát data. Stiskněte výběrové pole požadovaného datového formátu (příklad: MM/DD/RRRR zobrazí datum jako Duben 15, 2015) a stiskněte tlačítko Uložit.



Obraz 51: Nastavení formátu data

Čas: Pro nastavení času a formátu času stiskněte tlačítko Čas. Stiskněte výběrové pole 12hod. nebo 24hod., nastavte pomocí jednotlivých segmentů nastavovacího kolečka hodiny, minuty a AM/ FM (jen 12hod. formát) a stiskněte tlačítko Uložit.



Obraz 52: Nastavení času a zobrazení ve 12hod./24hod. formátu

Klidový režim: Stiskněte tlačítko Klidový režim, displej míchadla přejde po 15 minutách nečinnosti do klidového stavu. V klidovém režimu se na obrazovce zobrazuje výzva pro uživatele Aktivujte dotykem.



Obraz 53: Aktivace klidového režimu

Region: Pro nastavení regionu, ve kterém je zařízení provozováno, stiskněte tlačítko **Region**. Stiskněte textové pole Město, země a zahajte zadávání prvních písmen příslušného města. Po prvních třech znacích nabídne uživatelské rozhraní možné návrhy. Přijměte nabízený návrh, nebo pokračujte v zadávání celého názvu, pak stiskněte tlačítko **Uložit** na softwarové klávesnici.



Obraz 54: Nastavení regionu

Název zařízení: Pro pojmenování nebo přejmenování míchadla stiskněte tlačítko Název zařízení. Stiskněte textové pole Název zařízení a zahajte zadávání. Po dokončení stiskněte tlačítko Uložit na softwarové klávesnici.

< ♠	Sh 11 Ur	ake La :22:24/ hit Nam	b 1 AM e		(< ♠	Shake Lab 1 11:22:24AM Unit Name	Ø	
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	S	ave	Ē			
1	Z	Х	С	۷	В	Ν	М	@		1			_	
?!1	23							-	-				Save	0

Obraz 55: Pojmenování míchadla

Úprava menu: Stiskněte tlačítko Úprava menu, chcete-li upravit spodní dva symboly na hlavní navigační liště podle své potřeby. Přetáhněte požadovaný symbol z hlavní obrazovky na symbol, který chcete nahradit. Stiskněte tlačítko Uložit, změna se potvrdí.

UPOZORNĚNÍ Stisknutím tlačítka Reset menu můžete navigační lištu kdykoli resetovat do továrního nastavení.



Obraz 56: Úprava navigační lišty

Soubory a informace

Tato obrazovka obsahuje informace o sériovém čísle, verzi aplikačního softwaru grafického uživatelského rozhraní, verzi firmwaru hlavní řídicí jednotky míchadla, verzi souboru parametrů a informaci o aktuálním operačním systému.

Stisknutím tlačítka **Reset do továrního nastavení** můžete míchadlo uvést do továrního nastavení. Reset do továrního nastavení vyžaduje přístupové heslo administrátora, při resetu do továrního nastavení se smažou všechna nastavení, která byla provedena přes grafické uživatelské rozhraní. Protokol událostí se při resetu do továrního nastavení zachová (nesmaže).

< ☆ ■	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	<	Shake Lab 1 11:22:24AM Files and Info Main Controller 7000 0612 V3 Parameters 50154439 V1 Board Support Package: 4.1.15/jethro-15.0	î
Ē	Board Support Package: 4.1.15-jethro-15.0	Ē	Factory Reset	>
Ļ	Factory Reset	.	Firmware	\rightarrow
	(D		0

Obraz 57: Soubory a informace

Při rolování směrem dolů se zobrazí tlačítko **Firmware**, jak je znázorněno na pravé straně Obraz 57. Toto tlačítko můžete stisknout, chcete-li do míchadla nainstalovat nový firmware. Podrobné pokyny k instalaci nového firmwaru najdete v oddíle "4. 7. Instalace firmwaru" na straně 130.

UPOZORNĚNÍ Nový firmware je nutné získat od autorizovaného servisního technika.

<u>Servis</u>

Přístup k servisním nastavením je omezený a je dovolen jen autorizovaným servisním technikům.

3. 2. 4. Programy

Stiskněte symbol **Programy** na navigační liště, chcete-li zobrazit seznam programů nebo vytvořit nový program. Programy je možné vytvářet, upravovat, mazat, importovat a exportovat. Na Obraz 58 je znázorněna obrazovka Programy nového zařízení (vlevo) a již používaného zařízení (vpravo). V používaném zařízení je již k dispozici seznam uživatelem vytvořených programů.



Obraz 58: Programy

U každé položky seznamu programů je k dispozici symbol oka (), pomocí kterého otevřete náhled programu s podrobnostmi programu, a symbol tužky , pomocí kterého otevřete program pro editaci.

Nastavení preferencí pro všechny programy

Z obrazovky Programy přejdete tlačítkem **Rozšířená nastavení** na obrazovku **Rozšířená nastavení**. Na této obrazovce můžete určit, co se stane při spuštění programu. Tato nastavení jsou platná pro všechny programy, které vytvoříte.



Obraz 59: Nastavení preferencí programu

- 1. Stiskněte tlačítko **Rozšířená nastavení**.
- 2. Vyberte volbu pro Zobrazení časovače, kterou určíte, jak se má časovač programu zobrazovat:
 - » Uplynulý: jak dlouho už míchadlo běží od okamžiku, kdy jste stiskli tlačítko Start, nebo
 - » Zbývající: jak dlouho ještě míchadlo poběží, než uplyne čas nastavený na časovači.
- Chcete-li, aby se míchadlo rozběhlo ihned při spuštění programu, aktivujte volbu Spustit míchadlo při načtení programu. Standardně je pro spuštění programu nutné stisknout tlačítko Start.

UPOZORNĚNÍ Volba **Spustit míchadlo při načtení programu** funguje podle očekávání jen v případě, že je kryt míchadla zavřený. Ihned po zavření krytu se cyklus automaticky zahájí.

4. Stiskněte tlačítko Uložit.

Vytvoření programu

Je možné vytvořit a uložit až 99 programů.

- 1. Stiskněte tlačítko Vytvořit nový.
- 2. Zadejte název programu.



Obraz 60: Vytvoření programu

 Zadejte teplotu (jen míchadla s regulací teploty), otáčky a procesní čas programu v hodinách a minutách.

Pro deaktivaci regulace teploty u míchadla s regulací teploty nechte pole Teplota prázdné.

- Pro přidání dalšího kroku do programu rolujte obrazovkou směrem dolů a stiskněte tlačítko Přidat krok.
- 5. Stiskněte tlačítko Uložit, program se uloží.

Úprava programu

- 1. Stiskněte symbol tužky 🖍 vedle programu, který si přejete upravit.
- 2. Upravte požadovaná pole. Stiskněte tlačítko Uložit. Program se uloží s provedenými změnami.
- Můžete přidat další krok, a to stisknutím tlačítka Přidat krok na spodním okraji obrazovky. Jsou-li zobrazeny více než 3 kroky, rolujte obrazovkou směrem dolů, až uvidíte tlačítko Přidat krok.

UPOZORNĚNÍ Zobrazí-li se během míchacího procesu s více kroky alarmová nebo chybové hlášení, míchadlo se automaticky zastaví. Zobrazí-li se výstražné hlášení, míchací proces pokračuje.



Obraz 61: Přidat krok

Smazání programu

- 1. Stiskněte symbol tužky vedle programu, který si přejete upravit.
- 2. Stiskněte tlačítko Smazat. Zobrazí se okno, které vyzývá k potvrzení akce.



Obraz 62: Smazání programu

3. Stiskněte tlačítko **Smazat**. Zvolený program se smaže.

Provedení programu

- 1. Vyberte uložený program, který si přejete provést.
- 2. Chcete-li si program rychle prohlédnout, stiskněte symbol oka 💿 vedle programu.
- 3. Zobrazí se okno s rychlým náhledem programu.

Program Quick View		×	
Soluble Solution			
Step 1: 24°C / 300rpm , Step 2: 26°C / 350rpm , Step 3: 30°C / 500rpm ,	/ 1hr, 30min / 30min / 1hr		
Edit	Run		

Obraz 63: Rychlý náhled programu

4. Stiskněte tlačítko Spustit, program se načte do hlavní obrazovky.

V závislosti na zvoleném nastavení Automatické spuštění se program spustí okamžitě, nebo až po stisknutí tlačítka Start na hlavní obrazovce.

5. Po dokončení programu se zobrazí následující okno. Stiskněte tlačítko OK.



Obraz 64: Program dokončen

Import programů

Do míchadla je možné importovat programy vytvořené na jiném míchadle. Za tím účelem je do USB konektoru míchadla nutné připojit USB flash disk.

Následující obrazovka znázorňuje všechny programy nalezené na USB flash disku.



Obraz 65: Výběr programů pro import

Vyberte programy, které si přejete importovat. Stiskněte tlačítko Importovat.

Po úspěšném importu programů se zobrazí následující okno. Stiskněte tlačítko **OK**. Nyní můžete USB flash disk vysunout.



Obraz 66: Import dokončen

Export programů

Programy je možné exportovat z jednoho míchadla do druhého. Zkontrolujte, že je USB flash disk správně zasunutý.

Po zasunutí USB flash disku se zobrazí následující obrazovka:



Obraz 67: Výběr programů pro export

Vyberte programy, které si přejete exportovat. Stiskněte tlačítko Exportovat.

Po dokončení exportu se zobrazí následující okno. Stiskněte tlačítko **OK**. Nyní můžete USB flash disk vysunout.



Obraz 68: Export dokončen

Protokol událostí

Třetí symbol na navigační liště je Protokol událostí, který obsahuje datový blok s uživatelskými a systémovými událostmi. Obrazovka Protokol událostí se zobrazí po stisknutí symbolu Protokol událostí na navigační liště, jak je znázorněno na Obraz 69.



Obraz 69: Protokol událostí

Na této obrazovce se zobrazuje seznam aktuálních událostí, včetně data a času.

Výběrem jednotlivé události můžete zobrazit doplňující informace k dané události.

Události je možné filtrovat podle následujících kategorií:

- Alarmová hlášení
- Výstražná hlášení
- Změny nastavení
- Spuštění a zastavení
- Programové cykly
- Otevření krytu (jen míchadla s regulací teploty)

Po zvolení některého z filtrů se pravé tlačítko Zobrazit všechny změní na Filtr ZAP., jak je znázorněno na Obraz 70.

5/12/15	Unit Started		~
5/12/15 11:34:06AM	Unit Stopped		~
5/12/15 11:34:06AM	High Temp Alarm		~
5/12/15 11:34:06AM	Setpoint Change		~
Date/Time	Event	Filter ON	~
11:22:24AM Event Log			

Obraz 70: Filtrovaný protokol událostí

Export protokolu událostí

 Z rozbalovacího menu Exportované události vyberte události, které si přejete exportovat. Vyberte formát, do jakého se má protokol nebo zpráva exportovat.



Obraz 71: Výběr protokolu událostí pro export

2. Je možné zvolit předdefinovaný, nebo zákazníkem definovaný datový rozsah.

UPOZORNĚNÍ Exportovat je možné jen události z posledních šesti měsíců.

 Export se provádí pomocí USB flash disku. Aby bylo možné protokol nebo zprávu uložit, musí být zasunutý USB flash disk. Stiskněte tlačítko Exportovat protokol, protokol nebo zpráva se stáhne.



Obraz 72: Zasuňte USB flash disk pro export

<u>Grafy</u>

Na obrazovce Grafy se znázorňují údaje o otáčkách a teplotách (jen míchadla s regulací teploty) v čase. Na ose X se zobrazuje čas, na ose Y otáčky nebo teplota.

V teplotním grafu můžete zaznamenávat požadovanou, skutečnou a okolní teplotu v čase. Tlačítkem Upravit můžete zvolit, která ze 3 křivek se má do teplotního grafu zaznamenat. Dále můžete upravit zobrazovací rozsah na 1 den, 7 dní nebo vlastní sekvenci dní, a to buď celých 24 hodin, nebo jen ve vybraných hodinách každého dne. Tlačítkem **Aktualizovat** zaktualizujete grafické zobrazení podle provedených nastavení a zobrazíte všechny nové údaje, které se zaznamenaly od posledního načtení nebo poslední aktualizace grafu.

Showing:	5/12/15, 8AM 6	o 4PM	-	Se Se	tpoint	Actual	^	Showing: 5/1	2/15, 8AM to	4PM	Set	point	Actual	Ambie
350rpm			11					70.0°C						
300rpm								60.0°C						
250rpm							444	50.0°C						
200rpm	-							40.0°C						
150rpm								30.0°C		-			1 1	-
100rpm							F	20.0°C						
50rpm								10.0°C						
Orpm	8AM 5/12/15	9AM	10AM	11AM	12PM			0.0°C	8AM	9AM	10AM	11AM	1201	N

Obraz 73: Grafy otáček a teplot

Do USB konektoru míchadla můžete připojit USB flash disk, po stisknutí tlačítka **Exportovat** grafická data se grafické údaje o otáčkách a teplotách stáhnou.

Úprava grafu

1. Stiskněte tlačítko Upravit, chcete-li zahájit úpravu grafu.



Obraz 74: Úprava grafu: Výběr teploty nebo otáček pro záznam

- 2. Stisknutím tlačítka Teplota nebo Otáčky zvolte, který graf se má zobrazit.
- 3. Rolujte obrazovkou směrem dolů, až uvidíte volbu Datový rozsah.
- 4. Stiskněte 1 den, 7 dní nebo Vlastní, chcete-li vybrat vlastní datový rozsah.
- Zvolte pomocí jednotlivých segmentů nastavovacího kolečka datum zahájení a stiskněte tlačítko Od.

 Zvolte pomocí jednotlivých segmentů nastavovacího kolečka datum ukončení a stiskněte tlačítko Do.



Obraz 75: Výběr datového rozsahu grafu

- 7. Rolujte obrazovkou směrem dolů, až uvidíte volbu Časový rozsah.
- 8. Stiskněte 24hod. denně, nebo stiskněte Vlastní, chcete-li zobrazit určitý časový úsek a
 - Zvolte pomocí jednotlivých segmentů nastavovacího kolečka každodenní čas spuštění a stiskněte tlačítko Od.
 - Zvolte pomocí jednotlivých segmentů nastavovacího kolečka každodenní čas ukončení a stiskněte tlačítko Do.

UPOZORNĚNÍ Volbou časových bodů pro spuštění a ukončení se rozhodnete pro přesné zobrazení časového úseku mezi dvěma časovými body. Volbou 24 hod. denně se záznam zahájí každý den v 00:00 hod. a ukončí ve 23:59 hod.



Obraz 76: Výběr denních dob pro záznam

9. Stiskněte tlačítko Uložit, provedené změny nebo vlastní nastavení se uloží.

3.3. Příslušenství



POZOR

Zranění a biologické ohrožení rozbitými nádobami.

Neodborně nainstalované příslušenství může způsobit prasknutí skla a rozlití vzorků.

Zkontrolujte, že jsou díly příslušenství řádně nainstalované, použijte správné nářadí a šrouby.

Zkontrolujte, že díly příslušenství řádně sedí na základně.

Nádoby používejte vždy jen s příslušenstvím, které odpovídá jejich velikosti.



Nebezpečí pořezání o ostré hrany.

Se základnami a ostatním příslušenstvím zacházejte opatrně.

POZOR

Pohybující se základna může přiskřípnout prsty.

POZOR

UPOZORNĚNÍ Za řádnou instalaci příslušenství odpovídá sám zákazník.

Nesahejte prsty na/pod základnu v pohybu.

Není-li uvedeno jinak, platí popsané postupy instalace pro všechny varianty míchadel uvedené v tomto návodu.

Používejte vždy správné nářadí, především pak nářadí dodané s příslušenstvím. V případě ztráty některých z těchto položek můžete objednávat uvedené sady náhradních dílů ("Příslušenství" na straně 17). Používejte jen a výhradně uvedené nářadí a jen a výhradně dodané šrouby.

3. 3. 1. Montáž základny

POZOR Kompletní seznam kompatibilních základen pro každé míchadlo najdete v oddíle "1. 2. 1. Základny" na straně 18. Míchadla s regulací teploty je dovoleno provozovat jen se základnami předinstalovanými z výroby.

POZOR Nebezpečí pořezání o ostré hrany. Základnu uchopte při demontáži zespodu. Základnu nikdy nezvedejte za držáky baněk, které jsou k základně připevněné. Se základnami a ostatním příslušenstvím zacházejte opatrně.

POZOR V míchadle Solaris 4000 I / 4000 R hrozí nebezpečí přiskřípnutí a pohmoždění prstů a rukou při pohybující se základně 11x14. Pro míchadlo Solaris 4000 I / 4000 R nepoužívejte základnu 11x14.

POZOR Použijte výhradně šroubový materiál a upevňovací nářadí dodané se základnou. Použití jiných šroubů nebo nesprávného upevňovacího nářadí neumožňuje řádnou instalaci a může vést k poškození míchadla či příslušenství.

Základnu upevněte vždy všemi šrouby.

Použijte vždy správné upevňovací nářadí:

- 3/16" upevňovací nářadí (GT530066) pro všechny základny míchadel Solaris 2000, Solaris 2000 I/R a 4000 I/R
- 7/32" upevňovací nářadí (GT530080) pro všechny základny míchadel Solaris 4000

Univerzální základny

Každé míchadlo konstrukční řady Solaris 2000 a 4000 se dodává s univerzální základnou, šrouby a upevňovacím nářadím. Pro další aplikace je možné objednat další základny. Kompletní seznam kompatibilních základen pro každé míchadlo najdete v oddíle "1. 2. 1. Základny" na straně 18.

U míchadel 2000 I, 2000 R, 4000 I a 4000 R je už základna předinstalovaná v míchadle. Základna ani šrouby nejsou k dispozici jako samostatné díly.



Obraz 78: Univerzální základna, instalovaná na míchadle Solaris 2000 (vlevo) a míchadle Solaris 4000 (vpravo)

1. Opatrně položte základnu vodorovně na míchadlo a jeho 3 montážní body.

Základny pro míchadla Solaris 2000 mají 3 montážní otvory. Základny pro míchadla Solaris 4000 mají 4 montážní otvory.

Základny pro míchadla Solaris 2000 I/R a 4000 I/R mají 3 montážní otvory.

 Šrouby opatrně utáhněte, základna je připevněna k míchadlu. Utahování šroubu přerušte v okamžiku, kdy se upevňovací nářadí začne prohýbat.

Dvoupatrové základny

UPOZORNĚNÍ Dvoupatrové základny nejsou vhodné pro míchadla Solaris 2000 I/R a 4000 I/R.



Obraz 79: Konstrukční skupina dvoupatrové základny

Spojte horní a spodní základnu pomocí 4 vzpěr nainstalovaných ve všech rozích. Upevněte vzpěry příslušnými šrouby k horní a spodní základně.

Šrouby opatrně utáhněte, vzpěry jsou spojené se základnami. Utahování šroubu přerušte v okamžiku, kdy se upevňovací nářadí začne prohýbat.

Spodní základna disponuje montážními otvory pro připevnění k míchadlu. Zkontrolujte, že se spodní základna při sestavování nachází dole.

Příslušenství

UPOZORNĚNÍ Než na konstrukční skupinu dvoupatrové základny postavíte nádoby, proveďte poslední kontrolu a ujistěte se, že se konstrukční skupina nekývá.



Obraz 80: Dvoupatrová základna na míchadle Solaris 4000

 Opatrně položte smontovanou dvoupatrovou základnu vodorovně na míchadlo a jeho montážní desku.

Základny pro Solaris 2000 mají 3 montážní otvory, základny pro Solaris 4000 pak 4 montážní otvory.

 Šrouby opatrně utáhněte, základna je připevněna k míchadlu. Utahování šroubu přerušte v okamžiku, kdy se upevňovací nářadí začne prohýbat.

3. 3. 2. Nasazení držáků baněk a nádob

POZOR Biologické ohrožení rozbitými nebo netěsnými nádobami. Neodborně nainstalované příslušenství může vést k rozlití vzorků. Zkontrolujte, že jsou díly příslušenství řádně nainstalované, použijte správné nářadí a šrouby. Pro nasazení je nutné použít šroubovák (PH2) s délkou dříku 150 mm (6 palců) (art. č. 75004131). Zkontrolujte, že díly příslušenství řádně sedí na základně. Nádoby používejte vždy jen s příslušenstvím, které odpovídá jejich velikosti. Nádoby je nutné instalovat řádně a v neporušeném stavu.

POZOR Nebezpečí pořezání o ostré hrany. Se základnami a ostatním příslušenstvím zacházejte opatrně.

POZOR Bezpečné upevnění držáku baňky je zajištěné jen při použití dodaných šroubů. Vložte dodané šrouby do všech montážních otvorů držáku baněk.

UPOZORNĚNÍ Příliš dlouhé šrouby mohou ohrozit stabilitu a funkčnost vyměnitelné základny. Dbejte na to, aby byly pro připevnění použity jen dodané šrouby.

Držáky baněk

Každý držák baněk se skládá z přídržné spony, jedné nebo dvou pružin (podle velikosti držáku/ lahve) a šroubů pro připevnění k základně. Použijte výhradně šrouby dodané s držákem baněk.



① Pružinové pouzdro ② Pružina ③ Šrouby ④ Přídržná spona ⑤ Pružina

Obraz 81: Držák baněk se 2 pružinami

- 1. V případě potřeby připevněte na přídržné spony pružinu, jak je znázorněno na obrázcích.
- Pružinová pouzdra se nasazují mezi ramena přídržné spony, jak je znázorněno na obrázku. Některé přídržné spony používají dvě pružiny. Druhá pružina se nasazuje okolo spodní části přídržné spony, a to po připevnění držáku baněk k základně.
- 3. Připevněte přídržnou sponu k základně pomocí dodaných šroubů.

<u>Nádoba</u>

- Vsaďte požadovanou nádobu opatrně do držáku baněk. Nejprve dostatečně roztáhněte přídržnou pružinu, aby bylo dno nádoby možné umístit doprostřed držáku baněk. Poté nádobu zlehka zasuňte do správné pozice a upevněte ji ve spodní, širší části držáku baněk. Hrdlo nádoby drží v bezpečné pozici pružina.
- Před zapnutím zařízení zkontrolujte, že jsou všechny nádoby dobře upevněné.

Nádoba by měla být podle možností uzavřená zátkou, předejdete tak vystříknutí látek při míchání.

3. 3. 3. Instalace čtvercového držáku baněk

POZOR Biologické ohrožení rozbitými nebo netěsnými nádobami. Neodborně nainstalované příslušenství může vést k rozlití vzorků. Zkontrolujte, že jsou díly příslušenství řádně nainstalované, použijte správné nářadí a šrouby. Pro nasazení je nutné použít šroubovák (PH2) s délkou dříku 150 mm (6 palců) (art. č. 75004131). Zkontrolujte, že díly příslušenství řádně sedí na základně. Nádoby používejte vždy jen s příslušenstvím, které odpovídá jejich velikosti. Nádoby je nutné instalovat řádně a v neporušeném stavu.

POZOR Nebezpečí pořezání o ostré hrany. Se základnami a ostatním příslušenstvím zacházejte opatrně.



Čtvercový držák baněk

Každý čtvercový držák baněk se skládá ze základní desky s montážními otvory a čtyř pružinových přídržných spon. Součástí jsou také šrouby pro připevnění držáku baněk k základně. Použijte výhradně šrouby dodané s držákem baněk.



① Montážní otvory ② Pružina ③ Přídržná spona ④ Šrouby

Obraz 82: Čtvercový držák baněk

- 1. Zavěste případně všechny pružiny 2 do přídržných spon 3, jak je znázorněno na Obraz 82.
- 2. Položte skupinu s přídržnými sponami na základnu a slícujte montážní otvory ①.
- 3. Připevněte skupinu s přídržnými sponami k základně pomocí dodaných šroubů ④.
- 4. Vsaďte čtvercovou nádobu shora do držáku.
- 5. Zkontrolujte, že nádoba dosedá na základní desku a pevně sedí v držáku.

3. 3. 4. Instalace stojanu na zkumavky

POZOR Nebezpečí pořezání o ostré hrany. Se základnami a ostatním příslušenstvím zacházejte opatrně.

Každý stojan na zkumavky se skládá ze dvou montážních třmenů (drátěných nebo plechových) s jedním upevňovacím knoflíkem a dvěma šrouby s rýhovanou hlavou pro připevnění k základně. Každá konstrukční skupina stojanu na zkumavky se dodává v předmontovaném stavu vč. stojanu na zkumavky a pěnové vložky (pěnových vložek). Použijte vždy jen šrouby dodané se stojanem.



Obraz 83: Konstrukční skupina stojanu na zkumavky s montážními třmeny (drátěnými)



Obraz 84: Konstrukční skupina stojanu na zkumavky s montážními třmeny (plechovými)

- Nainstalujte montážní třmeny pomocí šroubů s rýhovanou hlavou na základnu. Utáhněte šrouby s rýhovanou hlavou pevně rukou.
- 2. Nasaďte stojan na zkumavky do montážních třmenů.
- Nastavte pomocí upevňovacích knoflíků požadovaný úhel. Utáhněte upevňovací knoflíky pevně rukou.

3. 3. 5. Instalace držáku stojanu na zkumavky s nastavitelným úhlem

POZOR Nebezpečí pořezání o ostré hrany. Se základnami a ostatním příslušenstvím zacházejte opatrně.

Držáky stojanů na zkumavky a stojany na zkumavky jsou k dispozici v různých velikostech. Seznam najdete v oddíle "1. 2. 5. Držáky stojanů na zkumavky s nastavitelným úhlem" na straně 28. Všechny držáky stojanů na zkumavky je možné nastavit do sedmi pozic, naklápět z 0° v obou směrech a aretovat v 15°, 30° a 45°.

Použijte vždy jen šrouby dodané s držákem stojanu.



Obraz 85: Držák stojanu na zkumavky s nastavitelným úhlem vč. nainstalovaného stojanu

- 1. Nainstalujte držák stojanu ① pomocí dodaných šroubů na základnu míchadla.
- Mírně rozepřete kovové spony ③ na obou koncích výklopné části ⑥ držáku a vsaďte do něj stojan na zkumavky ②.
- 3. Zajistěte stojan na zkumavky 2 opatrným uvolněním spon 3.
- Vytáhněte hlavy

 upevňovacích knoflíků na obou koncích držáku směrem ven. Knoflíky se odblokují otočením o 1/4 otáčky.
- 5. Natočte výklopnou část 6 držáku 1 do požadovaného úhlu 15°, 30° nebo 45°.
- Pro vyjmutí stojanu na zkumavky rozepřete kovové spony ③ na obou koncích výklopné části ⑥ držáku a vytáhněte stojan ① ven.

3. 3. 6. Instalace držáku pro mikrotitrační/DeepWell destičky

POZOR Nebezpečí pořezání o ostré hrany. Se základnami a ostatním příslušenstvím zacházejte opatrně.



DeepWell destička

② Šroub

③ Držák mikrotitrační destičky

Obraz 86: Konstrukční skupina držáku pro mikrotitrační/DeepWell destičky

- 1. Umístěte rám mikrotitrační destičky na základnu.
- Připevněte rám mikrotitrační destičky k základně pomocí šroubů dodaných se sadou mikrotitrační/DeepWell destičky.
- 3. Nasaďte mikrotitrační/DeepWell destičku do rámu.
- 4. Zkontrolujte opatrným přizvednutím pevný sed mikrotitrační/DeepWell destičky.

3. 3. 7. Instalace nastavitelného držáku nádob

POZOR Nebezpečí pořezání o ostré hrany. Se základnami a ostatním příslušenstvím zacházejte opatrně.



① Montážní otvory ② Stavitelné tyče ③ Upevňovací knoflíky ④ Boční stěna ⑤ Protiskluzová podložka
 ⑥ Šrouby

Obraz 87: Konstrukční skupina jednostupňového výškově nastavitelného držáku nádob

Příslušenství





① Montážní otvory ② Stavitelné tyče ③ Upevňovací knoflíky ④ Boční stěna ⑤ Protiskluzová podložka
 ⑥ Šrouby

Obraz 88: Konstrukční skupina dvoustupňového výškově nastavitelného držáku nádob

- 1. Nadzvedněte protiskluzovou podložku a odkryjte montážní otvory (poz. ① na Obraz 88).
- 2. Postavte nastavitelný držák na základnu a slícujte montážní otvory.
- 3. Připevněte nastavitelný držák k základně pomocí dodaných šroubů 6.
- 4. Položte protiskluzovou podložku (5) na dno nastavitelného držáku.
- 5. Na protiskluzovou podložku (5) postavte nádobu/nádoby.
- 6. Povolte upevňovací knoflíky 3.
- 7. Stavitelné tyče 2 umístěte tak, aby byly nádoby bezpečně upevněné.
- 8. Upevňovací knoflíky 3 opět utáhněte.
- 9. Opatrným přizvednutím zkontrolujte, že je nádoba bezpečně upevněná.

3. 3. 8. Instalace stojanu pro kádinky

POZOR Nebezpečí pořezání o ostré hrany. Se základnami a ostatním příslušenstvím zacházejte opatrně.



Obraz 89: Konstrukční skupina stojanu pro kádinky

- 1. Položte spodní desku 2 na základnu.
- 2. Na spodní desku položte stojan pro kádinky ①.
- 3. Slícujte montážní otvory se základnou.
- 4. Připevněte skupinu stojanu pro kádinky pomocí dodaných šroubů ④.

3. 3. 9. Instalace držáku pro dělicí nálevku

POZOR Nebezpečí pořezání o ostré hrany. Se základnami a ostatním příslušenstvím zacházejte opatrně.

UPOZORNĚNÍ Zkontrolujte, že používáte správné šrouby. Součástí dodávky jsou šrouby dvou typů. Delší šrouby s červeným zajišťovacím lakem jsou určené pro zajištění svislého stojanu. Kratší šrouby bez zajišťovacího laku jsou určené jen pro instalaci držáku pro dělicí nálevku na základnu.

Držák pro dělicí nálevku je na základnu míchadla možné instalovat ve dvou základních směrech: vodorovně a svisle.



Obraz 90: Instalační pozice držáku pro dělicí nálevku

Vodorovnou pozici ③ na Obraz 90 je možné výškově upravit a nálevku zpolohovat šikmo, jak je znázorněno na ④. Držák pro dělicí nálevku je dále možné zkombinovat se svislým stojanem (nabízeným jako příslušenství), který umožňuje nálevku zpolohovat svisle, jak je znázorněno na ① a ②. Držák je možné nainstalovat tak, aby kohout nálevky směřoval dolů ①, nebo obráceně ②.

Příprava svislé instalace



Obraz 91: Držák pro dělicí nálevku (vlevo) s příslušenstvím pro svislý stojan (vpravo)

- Upevněte držák pro dělicí nálevku na dlouhý konec svislého stojanu, jak znázorňují poz. ① a ② na Obraz 90.
- 2. Slícujte čtyři montážní otvory držáku nálevky s otvory svislého stojanu (2 na Obraz 91):
 - a. Pro instalaci kohoutu nálevky dole použijte otvory pro standardní svislou instalaci (3 na Obraz 91).
 - Pro instalaci kohoutu nálevky nahoře použijte otvory pro obrácenou svislou instalaci (@ na Obraz 91).
 - c. Upevněte držák pro dělicí nálevku pomocí dodaných šroubů (⑤ na Obraz 91) do svislého stojanu.

Montáž k základně

- 1. Položte konstrukční skupinu držáku pro dělicí nálevku na základnu.
- 2. Slícujte montážní otvory se základnou.
- Připevněte konstrukční skupinu držáku pro dělicí nálevku k základně pomocí šroubů dodaných s držákem (6 na Obraz 91 a Obraz 92).

Instalace nálevky



 Dělicí nálevka
 Zadní držák nálevky
 Přední držák nálevky
 Upevňovací knoflík, nastavení výšky
 Upevňovací knoflík, nastavení délky
 Šrouby bez zajišťovacího laku

Obraz 92: Konstrukční skupina držáku pro dělicí nálevku

- Povolte upevňovací knoflík pro nastavení délky (5) a posuňte přední držák nálevky (3) zcela dopředu.
- 2. Zaveďte zadní konec nálevky ① do zadního držáku nálevky ②.
- Posuňte přední držák nálevky ③ dozadu a vložte nálevkové hrdlo do otvoru předního držáku nálevky ③.
- 4. Nastavte přední držák nálevky ③ do vhodné výšky a zajistěte ho utažením upevňovacího knoflíku pro nastavení výšky ④. Úpravou výšky je dělicí nálevku možné zplohovat šikmo.
- 5. Zajistěte přední držák nálevky ③ utažením upevňovacího knoflíku pro nastavení délky ⑤.

3. 3. 10. Instalace univerzálního odkládacího tácu

POZOR Nebezpečí pořezání o ostré hrany. Se základnami a ostatním příslušenstvím zacházejte opatrně.



Obraz 93: Konstrukční skupina univerzálního odkládacího tácu

Při instalaci postupujte následujícím způsobem:

- 1. Položte univerzální odkládací tác 2 na základnu.
- 2. Slícujte montážní otvory se základnou.
- 3. Připevněte univerzální odkládací tác 2 pomocí šroubů dodaných s tácem 3.
- 4. Vložte do odkládacího tácu 2 gumovou podložku ①.
- 5. Položte na tác lahve a zajistěte je dodanými O-kroužky ④, viz Obraz 94.

UPOZORNĚNÍ O-kroužky doporučujeme používat pro otáčky nad 100 ot./min.



Obraz 94: Univerzální odkládací tác s upevněnými nádobami a O-kroužky

UPOZORNĚNÍ Dbejte na to, aby upevněné O-kroužky nebyly přepnuté. Při napínání O-kroužků dosáhnete v určitém okamžiku mrtvého bodu elasticity, za kterým se už O-kroužky nenapínají. Po dosažení tohoto bodu se O-kroužky nesnažte dále napínat, dojde k jejich přetržení.

3. 3. 11. Položení a použití adhezní podložky

Adhezní podložky jsou vhodné pro nádoby se širokým, plochým dnem, např. pro Erlenmeyerovy baňky, Fernbachovy baňky, Petriho misky, kultivační lahve, lahve na média, mikrotitrační destičky, odměrné baňky, kádinky. Obecně je možné používat všechny nádoby, které mohou stát zpříma. Se zmenšujícím se poměrem plochy dna vůči výšce se však snižují max. dosažitelné míchací otáčky. Vysoké nádoby s malým dnem jsou pro použití méně vhodné, viz Obraz 95. Velké Erlenmeyerovy baňky (např. 3000ml) přilnou lépe než malé (např. 100ml).



Obraz 95: Adhezní podložka: vhodné tvary baněk

Obecně nabízíme dva druhy adhezních podložek:

- Podložka s nízkou přilnavostí (transparentní) 75004111 a 75004117
 - » snadná manipulace díky nižší přilnavosti,
 - » vhodná především pro kádinky,
 - » vhodná pro všechny nádoby do max. otáček 200 ot./min (použití pro tkáňové kultury).
- Podložka s vysokou přilnavostí (černá, transparentní) 75004126 a 75004127
 - » spolehlivá fixace nádob díky vyšší přilnavosti,
 - » vhodná pro všechny nádoby,
 - » max. otáčky 400 ot./min (až do 500ml Erlenmeyer).

<u>Nářadí</u>

Aplikační váleček zobrazený na Obraz 96 se dodává jako příslušenství k adhezním podložkám. Slouží jako pomůcka pro přitlačení adhezní podložky k základně.



Obraz 96: Aplikační váleček pro adhezní podložky

Vložení suché podložky

- 1. Zkontrolujte, že je základna míchadla čistá a suchá.
- 2. Z jedné strany adhezní podložky odstraňte ochrannou fólii.
- 3. Nalepte podložku lepicí stranou směrem dolů na základnu.
- 4. Odstraňte ochrannou fólii z horní strany.

Vložení vlhké podložky

- 1. Z obou stran adhezní podložky odstraňte ochrannou fólii.
- 2. Namočte podložku do vody nebo ji důkladně opláchněte.
- 3. Nechte podložku okapat.
- 4. Rozprostřete vlhkou podložku na čistou základnu a umístěte ji podle potřeby.
- 5. Nechte podložku na základně 24 hodin uschnout, až pak podložku použijte.

Poznámky k vlastnostem nádob

Max. otáčky závisí na druhu nádoby. Nádoby se liší velikostí, tvarem, tvarem dna, materiálem, drsností povrchu a čistotou. Než začnete míchadlo používat, posuďte nejprve všechny uvedené parametry a zvolte nejvhodnější podložku.

Vhodnost laboratorního materiálu pro použití s adhezními podložkami shrnuje Tabulka 53.

Plastové / skleněné výrobky	Hodnocení přilnavosti	Kov	Hodnocení přilnavosti
Sklo (borosilikátové sklo)	5	Ocel, nerezová	5
Polypropylen (PP)	3	Hliník	5
Polypropylen kopolymer (PPCO)	3	Hliník eloxovaný	4
Polyethylen (HPDE) (LPDE)	3		
Polykarbonát (PC)	4		
Polyethylentereftalát (PET)	4		
Polyethylentereftalát	4		
kopolyester (PETG)			
Polymethylpenten (PMP)	3		
Polytetrafluorethylen (PTFE)	0*		
Tetrafluorethylen-	0*		
hexafluorpropylen kopolymer			
(FEP)			
Perfluoralkoxy polymer (PFA)	2		
Silikon	1*		
Hodnocení: 0 = žádná přilnavost; (PTFE, FEP a silikon).	5 = nejlepší přilnavost	t / * Nepoužívejte nádoby s	hodnocením 0 a 1

Tabulka 53: Kompatibilita materiálů nádob

UPOZORNĚNÍ Pamatujte, že max. míchacích otáček je možné dosáhnout jen s materiály, které jsou v Tabulka 53 hodnocené stupněm 5.

UPOZORNĚNÍ U jiných než skleněných materiálů proveďte vlastní zkoušky a určete, který druh materiálu je vhodný pro jaký náklad, otáčky, procesní čas a teplotu.
Rozsah otáček a mezní procesní čas pro skleněné nádoby (30% plnění)

Mezní hodnoty otáček a procesního času při teplotě okolí pro skleněné nádoby (30% plnění). Vyšší teploty snižují přilnavost. Při změně teploty se nádoby mohou uvolnit od podložky. Míchací proces spouštějte s aklimatizovanými nádobami.

	Mikrotitrační/DeepWell destičky	Kádinka					
	96 x 2 ml	25 ml	125 ml	250 ml	500 ml		
125 ot./min	neomezený procesní čas						
150 ot./min	72 h.	72 h.	72 h.	72 h.	72 h.		
200 ot./min	72 h.	72 h.	72 h.	72 h.	72 h.		
250 ot./min	72 h.						
300 ot./min							
350 ot./min							
400 ot./min							

Tabulka 54: Podložka s nízkou přilnavostí - rozsah otáček a mezní procesní čas při teplotě okolí

	Mikrotitrační/ DeepWell destičky	Erlenmeyer								
	96 x 2 ml	25 ml	125 ml	250 ml	500 ml	1000 ml	2000 ml	3000 ml	5000 ml	
125 ot./min	neomezený pro	omezený procesní čas								
150 ot./min	72 h.	72 h.	72 h.	72 h.	72 h.	72 h.	72 h.	72 h.	72 h.	
200 ot./min	72 h.	72 h.	72 h.	72 h.	72 h.	72 h.	72 h.	72 h.	72 h.	
250 ot./min	72 h.	72 h.	72 h.	72 h.	72 h.	72 h.	72 h.	48 h.	24 h.	
300 ot./min	72 h.	72 h.	72 h.	72 h.	72 h.	72 h.	48 h.	24 h.		
350 ot./min	48 h.	48 h.	48 h.	48 h.	48 h.	48 h.	24 h.			
400 ot./min	24 h.	24 h.	24 h.	24 h.	24 h.					

Tabulka 55: Podložka s vysokou přilnavostí - rozsah otáček a mezní procesní čas při teplotě okolí

<u>Příprava</u>

POZOR Při manipulaci se skleněnými výrobky noste vždy ochranné brýle a rukavice.

Před umístěním nádob na adhezní podložku:

- 1. Zkontrolujte nádoby, zda nejsou poškozené. Poškozené nádoby nepoužívejte.
- 2. Otřete nádobu čistou a suchou utěrkou, odstraňte veškerou vlhkost.
- 3. Odstraňte tuk a ostatní nečistoty.
- 4. Zkontrolujte, že jsou dno nádoby a adhezní podložka suché a čisté.

POZOR l jen sotva viditelný film nebo jediná kapka vody na dně nádoby může po určitém procesním čase způsobit ztrátu přilnavosti a uvolnění nádoby od podložky.



Obraz 97: Adhezní podložka: utření nádoby do sucha

5. Aby nedošlo k rozlití kapaliny, zmenšete plnicí množství nebo nádoby uzavřete.

POZOR I když nádoba k podložce bezvadně přilne, může jediná kapka kapaliny, která steče po vnější stěně nádoby, po určitém procesním čase způsobit uvolnění nádoby od podložky.

Vložení nádob

- 1. Umístěte nádoby na podložku a zatlačte je směrem dolů.
- 2. Opatrným tlakem/tahem zkontrolujte, že všechny nádoby k podložce spolehlivě přilnuly.

UPOZORNĚNÍ Pamatujte, že při nižších teplotách nebo delším provozu může docházet ke kondenzaci vody, v jejímž důsledku pak může dojít k neúmyslnému uvolnění nádob od podložky. Teplota vzorků a nádob by neměla překročit uvedený teplotní rozsah míchadla. Další informace k možným teplotním rozsahům míchadla najdete v oddíle "1. 1. Technické údaje" na straně 11.

Vyjmutí nádob

- 1. Stiskněte tlačítko Stop na displeji, nebo počkejte, než se cyklus dokončí.
- 2. Počkejte, než se základna zcela zastaví.
- Zatáhněte hrdlo nádoby opatrně a rovnoměrně do strany a počkejte několik sekund. U nádob s velkou plochou dna trvá uvolnění od podložky několik sekund.



Obraz 98: Adhezní podložka: opatrné odtlačení/odtažení nádoby

POZOR Nádobu tlačte/táhněte jemně a bez použití velké síly! Nádoby se mohou náhle uvolnit a může dojít k rozlití jejich obsahu!

4. Není-li nádobu možné uvolnit, nalijte okolo dna trochu vody.

Před dalším použitím nechte podložku vyschnout na vzduchu.

Odstranění podložky

 Uvolněte adhezní podložku nadzvednutím na jedné straně a pomalu ji stáhněte ze základny, jak je znázorněno na Obraz 99.



Obraz 99: Adhezní podložka: odstranění nadzvednutím a stažením

Čištění a dezinfekce

Při snížení přilnavosti podložku vyčistěte.

POZOR Nepoužívejte žádná jiná rozpouštědla, mycí ani čisticí prostředky než jemné mýdlo na nádobí bez přidaných prostředků pro péči o pokožku. Doporučená koncentrace je max. 2 ml prostředku na 1 litr vody.

Příslušenství

UPOZORNĚNÍ Přímý kontakt podložky s koncentrovaným čisticím prostředkem snižuje její přilnavost. V takovém případě není proplach dostatečným řešením. Po přímém kontaktu s čisticím prostředkem je podložku nutné propláchnout, namočit na 24 hodin do vody a pak nechat 24 hodin uschnout. Tímto způsobem je přilnavost možné plně obnovit.

UPOZORNĚNÍ Dezinfikujte jen 70% isopropanolem nebo 70% etanolem.

Lehké znečištění nebo zaprášení:

- 1. Podložku na základně otřete shora vlhkou utěrkou nebo podložku dezinfikujte.
- 2. Podložku navlhčete.
- 3. Setřete zbytky nečistot houbou.
- 4. Otřete podložku houbou navlhčenou v čisté vodě.
- 5. Podložku případně dezinfikujte.
- 6. Nechte podložku 24 hodin uschnout.

Silné znečištění:

- 1. Odstraňte podložku ze základny a kompletně ji vyčistěte.
- 2. Namočte podložku do vody se zředěným mýdlem na nádobí.
- 3. Odrhněte podložku důkladně houbou, odstraňte všechny zbytky nečistot.
- 4. Po vyčištění podložku důkladně propláchněte vodou.
- 5. Podložku případně dezinfikujte.
- 6. Nechte podložku 24 hodin uschnout.
- Vložte podložku podle popisu v oddíle "Vložení suché podložky" na straně 107 nebo "Vložení vlhké podložky" na straně 107.

3. 3. 12. Instalace rozdělovače plynu

Rozdělovač plynu umožňuje přivádět během míchacího provozu inertní plyn (např. dusík nebo oxid uhličitý) do více nádob současně (objednací informace najdete v oddíle "1. 2. 12. Obecné příslušenství" na straně 42).

Rozdělovač plynu je volitelné, továrně instalované příslušenství pro míchadla Solaris 2000 I/R a 4000 I/R. Není-li rozdělovač v zařízení nainstalovaný z výroby, je ho možné objednat a nainstalovat také dodatečně zákazníkem. Všechna míchadla typu Solaris 2000 I/R a 4000 I/R mají na boční stěně kabelovou průchodku (informace k jejímu umístění najdete v oddíle "1. 4. Přehled výrobku" na straně 44) pro protažení přívodní hadice laboratorního plynu ke konstrukční skupině rozdělovače. Rozdělovač plynu je vybaven jedním vstupním otvorem pro přívod plynu a osmi výstupními otvory pro nádoby.



① Přívodní hadice laboratorního plynu

② Konstrukční skupina rozdělovače plynu

3 Výstupní otvory s hadicemi

- ④ Vstupní otvor
- 5 Upevňovací šrouby

Obraz 100: Rozdělovač plynu instalovaný v míchadle Solaris 4000 R

Při instalaci postupujte následujícím způsobem:

- 1. Připevněte rozdělovač plynu 2 k zadní stěně míchací komory pomocí dodaných šroubů 5.
- Odstraňte plastovou zátku z vnější strany kabelové průchodky (informace k jejímu umístění najdete v oddíle "1. 4. Přehled výrobku" na straně 44).
- 3. Do membrány gumové průchodky uvnitř míchací komory vyřízněte pomocí lámacího nože kříž.
- 4. Přiřízněte kus hadice na vhodnou délku a použijte ji jako přívodní hadici laboratorního plynu ①.
- 5. Upevněte přívodní hadici laboratorního plynu ① do háku vstupního otvoru ④ rozdělovače.
- Dbejte na to, aby se přívodní hadice laboratorního plynu ① nedostala do kontaktu s pákou krytu a protáhněte ji gumovou průchodkou směrem ven.

Příslušenství

- 7. Připojte přívodní hadici laboratorního plynu ① k přívodu plynu.
- Uřízněte osm hadicových kusů a připevněte je k výstupům ③ rozdělovače. Je-li nádob víc než osm, použijte Y-spojky; je-li nádob méně než osm, zasvorkujte nepoužívané hadice tak, aby nedocházelo ke ztrátě plynu.
- 9. Pro zavedení hadic do nádob použijte vhodné zátky a těsnění.

3. 4. Nakládání a použití v souladu s určením



Nebezpečí požáru v důsledku vyvolaných chemických reakcí.

Neuvádějte míchadlo do provozu s otáčkami, při kterých dochází k vystříknutí obsahu nádob.

VÝSTRAHA

Otáčky zvyšujte pomalu. Před použitím chemikálií doporučujeme provést zkušební provoz s vodou.



Neodborné naložení a poškozené příslušenství mohou ohrozit bezpečnost.

POZOR

- Zkontrolujte, že je náklad (příslušenství a vzorky) rozmístěný symetricky ke středu základny. Při použití dvoupatrové základny dbejte na to, aby byla spodní základna naložená víc než základna horní.
- Míchadlo nikdy nepřetěžujte. Maximální náklad viz "Technické údaje" na straně 11. Náklad zahrnuje hmotnost základny instalované na míchadle, příslušenství a vzorků.
- Před uvedením míchadla do provozu zkontrolujte, že je příslušenství řádně nainstalované. Postupujte podle pokynů v oddíle "Příslušenství" na straně 88.



POZOR

Nebezpečí zranění prstů nebo rukou při pádu/přiklopení krytu.

 Kryt míchadla se může neúmyslně přiklopit a může přivřít nebo přiskřípnout prsty nebo ruce a způsobit jejich zranění. K tomu může dojít v případě, že je pružina krytu vadná a není s to udržet kryt ve zcela otevřené pozici. Neprovozujte míchadlo s vadnou pružinou krytu. Kontaktujte zákaznický servis společnosti Thermo Fisher Scientific a nechte pružinu krytu ihned vyměnit.

Vkládání

Zkontrolujte, že je náklad (příslušenství a vzorky) rozmístěný symetricky ke středu základny. Při použití dvoupatrové základny bez využití celé její kapacity: pro dosažení optimálních výsledků naložte přednostně spodní základnu a náklad rozmístěte symetricky ke středu základny.

Při použití míchadla s regulací teploty musí být náklad rozmístěný tak, aby byl při zavírání krytu i při zavřeném krytu k dispozici dostatek místa, aby nemohlo dojít ke kolizi a poškození nákladu.

Obraz 101: Příklad správně naložené základny

												л÷ч,		
(• •					*	*	*	÷	*	- f			
1	• •			*		*	٠	*	٠	٠		÷	* (•))) 👻
ŀ	• •	•	*	*	٠	*	*	*	*	*	*	٠	. 🖷	
	• •		*	*	*	*	*	*	٠	*	٠	٠	· P	⊾{(•)}
ŀ	• •				*	*	*			*	•	*	H. •)} ❤
ŀ	• •	•	•	*	*	*	*	*	٠	*	٠	٠	, H	
ŀ					*		*	*	*	*	*	*	Ē	s ((•))
ŀ	• •		•	*		*	٠	*		š	٠		- (-))) 👻
1	• •	•		*		*	*	*	*	*	*	e	. 9	
ŀ	•			•		٠	*	*	*	*	٠		· P	⊾ {(•))
Ŀ	• •		•			*					*	e	*(*)) -

Obraz 102: Příklad chybně naložené základny

UPOZORNĚNÍ Maximální náklad zahrnuje hmotnost základny instalované na míchadle, příslušenství a vzorků. Další informace k maximálnímu nákladu najdete v oddíle "1. 1. Technické údaje" na straně 11.

Použití v souladu s určením

Informace k mezním hodnotám otáček/nákladu při použití výrobku slouží jen jako orientační. Klidný provoz s malými nebo žádnými vibracemi závisí na mnoha faktorech, např. na plnicím množství, typu a stavu držáků baněk a podložek, typu nádob, uspořádání nádob na základně, podkladu míchadla a požadovaných otáčkách.

Při vytváření nové nebo neznámé kombinace, nebo při změně některého z parametrů vždy platí: Otáčky zvyšujte pomalu nad 175 ot./min a kontrolujte, zda nedochází k nežádoucím vibracím nákladu ani pohybům míchadla. Za bezpečný provoz míchadla odpovídá sám zákazník.

Míchadla typové řady Solaris jsou speciálně konstruovaná pro nízký výdej tepla a nízké vibrace, a jsou proto velmi vhodná pro použití v celé řadě laboratorních zařízení, vč. klimatizačních komor, inkubátorů a laboratorních chladicích skříní. Vzhledem k různým/specifickým podmínkám okolí, např. stabilitě základen a držáků, nese odpovědnost za bezpečný provoz při každém použití v laboratorních zařízeních sám uživatel. Otáčky zvyšujte pomalu a věnujte pozornost neobvyklým pohybům zařízení.

Bezpečné otáčky

Bezpečné otáčky míchadel konstrukční řady Solaris v závislosti na nákladu najdete v Tabulka 56. Náklad je souhrn hmotností základny instalované na míchadle, příslušenství a vzorků. Další informace k maximálnímu nákladu najdete v oddíle "Technické údaje" na straně 11.

Тур	Max. bezpečné otáčky
Solaris 2000	175 ot./min
Solaris 4000	175 ot./min
Solaris 2000 I	250 ot./min
Solaris 2000 R	250 ot./min
Solaris 4000 I	250 ot./min
Solaris 4000 R	250 ot./min

Tabulka 56: Max. bezpečné otáčky

Maximální otáčky

Při max. otáčkách 525 ot./min je dovoleno provozovat míchadla konstrukční řady Solaris s omezeným nákladem podle Tabulka 57. Náklad je souhrn hmotností základny instalované na míchadle, příslušenství a vzorků. Zohledněte přípustné rozsahy otáček použitých držáků a nádob, které s těmito otáčkami případně nejsou kompatibilní.

Тур	Max. bezpečný náklad
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

Tabulka 57: Maximální náklad při maximálních otáčkách

UPOZORNĚNÍ Vždy pamatujte na to, že se držáky s rostoucími otáčkami začínají otevírat a vzhledem k intenzivnějšímu pohybu nádob způsobují další dočasné vibrace systému.

Grafy pro hmotnost/otáčky

Následující grafy pro hmotnost/otáčky vám mají pomoci definovat optimální kombinace otáček/ nákladu pro použití výrobku v souladu s určením. Zelená oblast znázorňuje kombinace otáček/ nákladu, které při míchání způsobují malé nebo žádné vibrace systému. Červená oblast znázorňuje kombinace otáček/nákladu, které při míchání způsobují silné vibrace a které mohou vést k nežádoucímu pohybu míchadla. Protože jsou tyto grafy jen orientační, je nutné dávat pozor především v těch fázích, kdy se zvolená kombinace otáček/nákladu blíží červené oblasti.

Pamatujte, že vaše specifické podmínky použití mohou způsobit nežádoucí chování nákladu či zařízení před dosažením znázorněné hraniční linie. Otáčky zvyšujte v krocích, aby bylo možné sledovat chování specifického nákladu. Náklad je souhrn hmotností základny instalované na míchadle, příslušenství a vzorků.

UPOZORNĚNÍ Vodorovné modré čáry znázorňují v následujících grafech hmotnost příslušné prázdné základny.



Solaris 2000

Obraz 103: Solaris 2000 – použití v souladu s určením

Solaris 4000



Obraz 104: Solaris 4000 – použití v souladu s určením



Solaris 2000 I / 2000 R

Obraz 105: Solaris 2000 I / 2000 R – Použití v souladu s určením

Nakládání a použití v souladu s určením



Solaris 4000 I / 4000 R

Obraz 106: Solaris 4000 I / 4000 R – Použití v souladu s určením

4. Údržba a péče



VÝSTRAHA

Nebezpečí při zacházení s nebezpečnými látkami.

- Při míchání nebezpečných materiálů dodržujte doporučení Příručky o biologické bezpečnosti v laboratořích (Laboratory Biosafety Manual zpracovaný Světovou zdravotnickou organizací WHO) a další platná národní doporučení. Při míchání mikrobiologických vzorků rizikové skupiny II (dle Příručky o biologické bezpečnosti v laboratořích zpracované Světovou zdravotnickou organizací WHO) je nutné používat aerosoltěsná biotěsnění. Příručku o biologické bezpečnosti v laboratořích Laboratory Biosafety Manual najdete na internetové stránce Světové zdravotnické organizace (www.who.int). U materiálů vyšší rizikové skupiny je nutné provést další bezpečnostní opatření.
- Po kontaminaci míchadla nebo dílů toxiny nebo patogenními látkami je nutné provést vhodná dezinfekční resp. dekontaminační opatření ("Dekontaminace" na straně 124, "Dezinfekce" na straně 124).
- Nastane-li nebezpečná situace, přerušte napájení míchadla a ihned opusťte prostor.



VÝSTRAHA

Poškození zdraví infekčními látkami.

Proniknou-li nedopatřením kapaliny nebo jiné materiály pod základnu, ihned míchadlo vypněte, vytáhněte zástrčku z elektrické sítě a sejměte základnu ("Základny" na straně 18).

Utřete rozlitou kapalinu v souladu se standardním laboratorním postupem. Použijte vhodné osobní ochranné prostředky.

Všechna interní nastavení a opravy musí provádět autorizovaný servisní technik. Uživatel nesmí otevírat plášť míchadla.

Dodržujte všechny pokyny k řádnému provádění údržby a čištění, které najdete v produktových informacích dodaných s příslušenstvím. Informace uvedené v následujícím oddíle slouží jen jako vodítko.

4.1. Základní principy

V zájmu zajištění ochrany osob, životního prostředí a materiálu jste povinni míchadlo a jeho příslušenství pravidelně čistit a v případě nutnosti dezinfikovat.

Společnost Thermo Fisher Scientific doporučuje provést vyčištění a ruční dezinfekci míchadla alespoň jednou měsíčně. Okolní vzduch obsahuje tisíce cirkulujících mikroorganizmů, které se usazují v míchadle a které mohou ohrozit zpracovávané kultury.

Společnost Thermo Fisher Scientific doporučuje používat 70% etanol, 70% isopropanol nebo dezinfekční prostředky na bázi kvartérních amoniových solí (10% nebo méně).

UPOZORNĚNÍ Mechanizmus mohou poškodit pronikající kapaliny. Zabraňte proniknutí kapalin, především organických rozpouštědel, do mechanizmu nebo uložení mechanizmu. Organická rozpouštědla rozkládají tuk v uložení mechanizmu. Proniknou-li kapaliny do částí míchadla, které nejsou pro vyčištění přístupné, neuvádějte míchadlo do provozu a kontaktujte zákaznický servis.

UPOZORNĚNÍ Neschválené postupy nebo prostředky mohou napadat materiály míchadla a způsobit chybnou funkci. Nepoužívejte jiné než zde popsané čisticí a dekontaminační postupy, jestliže si nejste jisti, že je alternativní postup pro dané materiály vhodný. Používejte jen čisticí prostředky, které nepoškozují použité materiály. V případě pochybností kontaktujte výrobce čisticího prostředku.

UPOZORNĚNÍ U míchadel s regulací teploty ponechte otevřený kryt, aby se mohl čisticí prostředek a voda odpařit.

- Před zahájením čištění, dezinfekce nebo dekontaminace vytáhněte zástrčku z elektrické sítě.
- Před zahájením čištění, dezinfekce nebo dekontaminace demontujte z míchadla nainstalované příslušenství a základnu (základny).
- K čištění materiálů použijte teplou vodu s jemným čisticím prostředkem a měkkou utěrku.
 V případě pochybností kontaktujte společnost Thermo Fisher Scientific. Nakonec opláchněte čistou vodou a důkladně vysušte.
- Nikdy nepoužívejte žíravé čisticí prostředky jako kyselinu fosforečnou, bělicí louh nebo čisticí prášky.
- Používejte jen dezinfekční prostředky s hodnotou pH 6–8.
- Rozlité kapaliny ihned odstraňte utěrkou nepouštějící vlákna, navlhčenou v nekorozivním čisticím prostředku (viz pokyny výrobce čisticího prostředku).
- Rozlitá kapalina může zatéct pod základnu. Pronikne-li rozlitý obsah pod základnu, demontujte základnu a důkladně vyčistěte misku pod základnou. Proveďte případně dezinfekci nebo dekontaminaci.
- Zkontrolujte komponenty míchadla a odstraňte rozlitá/usazená média nebo nečistoty.

Kontrola příslušenství

UPOZORNĚNÍ Nepoužívejte míchadlo ani příslušenství vykazující známky poškození. Pro zajištění bezpečnosti doporučujeme příslušenství nechat kontrolovat v rámci pravidelné údržby.

Po důkladném vyčištění příslušenství také zkontrolujte, zda není poškozené, opotřebené nebo zkorodované.

Kovové díly

Vykazuje-li příslušenství známky poškození, jako je koroze, opotřebení nebo praskliny, ihned ho vyřaďte z provozu.

Plastové díly

Zkontrolujte tyto díly, zda nevykazují známky blednutí, praskliny, trhliny nebo poškrábání. Objevíte-li známky poškození, vyřaďte zkoumaný díl ihned z provozu.

Adhezní podložky

Adhezní podložky zkontrolujte před každým použitím, zda nevykazují známky opotřebení nebo poškození, jako jsou praskliny, trhliny, zabarvení, žloutnutí, zkřehnutí, deformace, povrchové odření nebo chemické poškození. Adhezní podložky, které vykazují známky opotřebení nebo poškození, ihned zlikvidujte.

4.2. Čištění

UPOZORNĚNÍ Před zahájením čištění ověřte u výrobce čisticího prostředku, že čisticí prostředek nepoškozuje použité materiály.

UPOZORNĚNÍ Mechanizmus mohou poškodit pronikající kapaliny. Zabraňte proniknutí kapalin, především organických rozpouštědel, do mechanizmu nebo do uložení mechanizmu. Organická rozpouštědla rozkládají tuk v uložení mechanizmu. Vnější strany zařízení umyjte jemným mýdlovým roztokem a měkkou utěrkou, opláchněte čistou vodou a důkladně vysušte.

Další informace k řádnému čištění míchadla a použitého příslušenství najdete v oddíle "Základní principy" na straně 122.

UPOZORNĚNÍ U míchadel s regulací teploty ponechte otevřený kryt, aby se mohl čisticí prostředek a voda odpařit.

Ovládací panel

Ovládací panel vyčistěte následujícím způsobem:

- 1. Vytáhněte zástrčku z elektrické sítě.
- 2. Vyčistěte ovládací panel suchou utěrkou z mikrovláken.
- 3. V případě nutnosti utěrku z mikrovláken navlhčete vodou a ovládací panel znovu vyčistěte.
- 4. Nechte všechny komponenty vyschnout na vzduchu.

4.3. Dezinfekce

VÝSTRAHA Nebezpečí při zacházení s nebezpečnými látkami. Nedotýkejte se infikovaných dílů. Při kontaktu s kontaminovanými díly může dojít k těžké infekci. Infekční materiál může při rozbití nádoby nebo rozlití proniknout do míchadla. V případě kontaminace zajistěte, aby nebyly ohroženy další osoby. Zasažené díly ihned vydezinfikujte.

UPOZORNĚNÍ Použitím nevhodných dezinfekčních metod nebo prostředků může dojít k poškození materiálů. Zkontrolujte, že dezinfekční metoda nebo prostředek nepoškozuje použité materiály. V případě pochybností kontaktujte výrobce dezinfekčního prostředku. Dodržte bezpečnostní pokyny a pokyny k použití aplikovaných dezinfekčních prostředků.

- Otřete všechny komponenty a plochy 70% etanolem, podle potřebné intenzity dezinfekce. Přístupné elektronické komponenty nesmí přijít do kontaktu s kapalinou.
- 2. Nechte všechny komponenty vyschnout na vzduchu.

UPOZORNĚNÍ U míchadel s regulací teploty ponechte otevřený kryt, aby se mohl čisticí prostředek a voda odpařit.

Uživatel sám odpovídá za to, že stupeň dekontaminace splňuje jím stanovené požadavky.

4.4. Dekontaminace

VÝSTRAHA Nebezpečí při zacházení s nebezpečnými látkami. Nedotýkejte se kontaminovaných dílů. Při kontaktu s kontaminovanými díly může dojít ke kontaminaci. Kontaminovaný materiál může při rozbití nádoby nebo rozlití proniknout do míchadla. V případě kontaminace zajistěte, aby nebyly ohroženy další osoby. Zasažené díly ihned dekontaminujte.

UPOZORNĚNÍ Použitím nevhodných dekontaminačních metod nebo prostředků může dojít k poškození materiálů. Zkontrolujte, že dekontaminační metoda nebo prostředek nepoškozuje použité materiály. V případě pochybností kontaktujte výrobce dekontaminačního prostředku. Dodržte bezpečnostní pokyny a pokyny k použití použitých dekontaminačních prostředků.

Společnost Thermo Fisher Scientific doporučuje následující metodu.

- Otřete všechny komponenty a plochy 70% etanolem. Přístupné elektronické komponenty nesmí přijít do kontaktu s kapalinou.
- 2. Nechte všechny komponenty vyschnout na vzduchu.

UPOZORNĚNÍ U míchadel s regulací teploty ponechte otevřený kryt, aby se mohl čisticí prostředek a voda odpařit.

Uživatel sám odpovídá za to, že stupeň dekontaminace splňuje jím stanovené požadavky.

4.5. Autoklávování

UPOZORNĚNÍ Nikdy nepřekročte přípustné hodnoty pro teplotu a délku autoklávování. Použití chemických přísad v páře je zakázáno.

Před zahájením autoklávování:

- Demontujte všechny díly příslušenství.
- Důkladně opláchněte všechny stopy chemikálií a čisticích prostředků. Proveďte závěrečný oplach destilovanou vodou.

Informace k autoklávování najdete v následující tabulce:

Artikl	Možnost autoklávování	Technické údaje
Míchadlo	Ne	-
Základna	Ano	121 °C, 20 min
Držák baněk (bez pružinových pouzder)	Ano	121 °C, 20 min
Šroubovák	Ne	-
Držáky baněk pro mikrotitrační/DeepWell destičky	Ano	121 °C, 20 min
Adhezní podložka	Ne	-
Stojan pro kádinky	Ano	121 °C, 20 min
Držáky pro dělicí nálevky	Ano	121 °C, 20 min
Stojan na zkumavky – kovové díly	Ano	121 °C, 20 min
Stojan na zkumavky – plastové díly	Ne	-
Stojan na zkumavky – pěnové vložky	Ne	-
Nastavitelný držák nádob – kovové díly	Ano	121 °C, 20 min
Nastavitelný držák nádob – protiskluzová podložka	Ne	-
Univerzální odkládací tác – kovový tác	Ano	121 °C, 20 min
Univerzální odkládací tác – gumová podložka	Ne	-
Držák stojanu na zkumavky s nastavitelným úhlem – kovové díly	Ano	121 °C, 20 min
Držák stojanu na zkumavky s nastavitelným úhlem – plastové díly	Ne	-
Držák stojanu na zkumavky s nastavitelným úhlem – pěnové vložky	Ne	-

Tabulka 58: Možnost autoklávování jednotlivých materiálů

Zajistěte, aby byla zajištěna potřebná sterilita v souladu s interními požadavky.

Po autoklávování může dojít k mírné změně vzhledu a barvy.

4. 6. Kalibrace teploty

Požadovanou teplotu míchadla s regulací teploty je možné zkalibrovat pomocí teplotního měřicího přístroje s potřebnou přesností měření. Takto zkalibrovaná hodnota zaručuje úplnou reprodukovatelnost inkubovaného (jen vyhřívaného) nebo chlazeného (chlazeného a vyhřívaného) míchacího procesu.

4. 6. 1. Osvědčené postupy kalibrace teploty

Před kalibrací a během kalibrace teploty míchadla postupujte přesně podle uvedených pokynů. Každá chyba nebo odchylka od doporučeného postupu může vést k tomu, že se teplota vzorku nestabilizuje a kalibrační proces se nezdaří.

Osvědčené postupy jsou:

- Použijte baňku s objemem 250 ml.
- Naplňte baňku přesně 100 ml vody.
- Zavřete úplně kryt a nechte ho zavřený po celou dobu kalibračního procesu.
- Nechte uplynout dostatek času, aby bylo jisté, že se teplota vzorku zcela stabilizovala.
- Zajistěte, aby se podmínky okolí (teplota a vlhkost) během kalibračního procesu neměnily.
 I malé změny, jako otevření dveří/okna nebo změna nastavení klimatizace, mohou negativně ovlivnit výsledek probíhajícího kalibračního procesu.
- Kalibraci míchadla doporučujeme provádět jednou ročně.

4. 6. 2. Provedení kalibrace teploty

Pro provedení kalibrace postupujte následujícím způsobem:

- 1. Kalibraci je možné zahájit třemi způsoby:
 - » Na konci procesu základního nastavení popsaného v oddíle "Základní nastavení" na straně 54 můžete místo ukončení procesu provést kalibraci teploty.
 - » Míchadlo zobrazí připomenutí roční kalibrace.
 - » Kalibraci pak můžete zahájit stisknutím symbolu Nastavení na navigační liště a pak tlačítka Kalibrace dole na zobrazené obrazovce Kalibrace teploty.
 - » Kalibraci můžete zahájit podle potřeby stisknutím symbolu Nastavení na navigační liště a pak tlačítka Kalibrace.

UPOZORNĚNÍ Připomenutí kalibrace je možné deaktivovat, a to označením zaškrtávacího pole Deaktivace kalibračních hlášení na straně Nastavení. 2. Počkejte, než se na obrazovce Kalibrace teploty zobrazí startovací obrazovka, viz Obraz 107.



Obraz 107: Kalibrace: startovací obrazovka

3. Stiskněte tlačítko Start kalibrace, proces se zahájí.



Obraz 108: Kalibrace: výzva k naplnění Erlenmeyerovy baňky

- 4. Postupujte podle pokynů na obrazovce: Naplňte 250 ml Erlenmeyerovu baňku 100 ml vody.
- 5. Vložte baňku do držáku baněk a upevněte držák přibližně doprostřed základny míchadla.
- Zavěste do baňky teplotní snímač měřicího přístroje s potřebnou přesností měření. Snímač by měl být ponořený do kapaliny, neměl by se však dotýkat dna ani stěn baňky.
- 7. Uzavřete baňku, aby nedošlo k rozlití kapaliny.
- 8. Zavřete kryt nad základnou míchadla.

9. Stiskněte tlačítko **Pokračovat**.



Obraz 109: Kalibrace: výzva k zahájení kalibrace

UPOZORNĚNÍ Míchadlo zahájí míchací proces, jakmile stisknete tlačítko Start.

 Postupujte podle pokynů na obrazovce: Stiskněte tlačítko Start, chcete-li použít přednastavenou teplotu, nebo stiskněte pole Pož. teplota a zadejte jinou požadovanou hodnotu, až pak stiskněte tlačítko Start.

Na obrazovce se zobrazuje, jak se míchací komora ohřívá (nebo ochlazuje) na zvolenou kalibrační teplotu.



Obraz 110: Kalibrace: stabilizace komory na kalibrační teplotu

11. Počkejte alespoň dvě a půl hodiny (150 minut), než teplota v komoře dosáhne stabilního stavu.

Stabilní stav oznámí hlášení **Teplota v komoře se stabilizovala**, jak je znázorněno na Obraz 111 dole.



Obraz 111: Kalibrace: úspěšná stabilizace kalibrační teploty v komoře

- 12. Postupujte podle pokynů na obrazovce:
 - » Jakmile se zobrazí obrazovka na Obraz 111, pravidelně porovnejte zobrazenou teplotu s teplotou odečtenou teplotním snímačem. Považujete-li teplotu za stabilní, můžete stisknout tlačítko Upravit pro její korekci, nebo tlačítko Dokončit pro ukončení kalibračního procesu.



Obraz 112: Kalibrace: korekce teploty zobrazené míchadlem

- Pomocí zobrazeného nastavovacího kolečka proveďte korekci teploty zobrazené na míchadle na hodnotu zjištěnou měřicím přístrojem.
- 14. Stiskněte tlačítko Uložit, provedené změny se uloží.

15. Hlášení Kalibrace teploty je dokončena informuje o tom, že byl proces úspěšně dokončen.

<	Shake Lab 1 11:22:24AM Temp Calibration	
¢.	Step 5 of 5	
	Temperature calibration complete.	
	Re-Calibrate	G

Obraz 113: Kalibrace: úspěšné dokončení kalibrace teploty

Stisknutím tlačítka Znovu zkalibrovat můžete kalibrační proces zopakovat.

4. 6. 3. Vyhledání kalibrační teploty

Aktuální kalibrační teplotu je možné kdykoli vyhledat:

- Na navigační liště stiskněte symbol Nastavení a pak tlačítko Kalibrace dole na zobrazené obrazovce Kalibrace teploty.
- Zobrazí se startovací obrazovka kalibrace, viz Obraz 107. Na této obrazovce je uvedena aktuální Kalibrační teplota míchadla.

UPOZORNĚNÍ Alternativně můžete u míchadel s regulací teploty najít kalibrační teplotu na obrazovce, na které se nastavuje teplota. Příklad viz oddíl "Nastavení teploty" na straně 64.

4.7. Instalace firmwaru

Stisknete-li na obrazovce Soubory a informace (viz "Soubory a informace" na straně 77) tlačítko Firmware, zobrazí se výzva k zadání přístupového hesla administrátora a pak obrazovka Firmware znázorněná na Obraz 114. Na této obrazovce můžete provést instalaci nového firmwaru z USB flash disku.

V závislosti na uzavřené servisní smlouvě obdržíte od autorizovaného technika odkaz pro stažení nového firmwaru. Vždy instalujte jen firmware z autorizovaného zdroje. Použitím softwaru z neautorizovaných zdrojů může zaniknout záruka poskytnutá společností Thermo Fisher Scientific.

- Z odkazu získaného od autorizovaného technika si stáhněte komprimovaný soubor ve formátu ZIP.
- 2. Rozbalte ZIP soubor na prázdný USB flash disk.

UPOZORNĚNÍ Zkontrolujte, že je na USB flash disku dostatek místa pro rozbalení ZIP souboru.

- 3. Jestliže míchadlo ještě běží, zastavte ho stisknutím tlačítka Stop na displeji.
- Potvrďte všechna případná alarmová hlášení, jak je popsáno v oddíle "Provozní stav" na straně 66.

UPOZORNĚNÍ Pokusíte-li se provést aktualizaci při běžícím míchadle, zobrazí se následující hlášení: **Aktualizaci není možné provést, jestliže se míchadlo nachází v míchacím procesu.** Součástí tohoto hlášení je tlačítko **Zastavit míchadlo**, pomocí kterého můžete proces ukončit, pak můžete provést aktualizaci.

- 5. Na navigační liště stiskněte symbol Soubory a informace.
- 6. Rolujte obrazovkou Soubory a informace až dolů, zobrazí se tlačítko Firmware.
- Stiskněte tlačítko Firmware, otevře se obrazovka Firmware, na které najdete aktuálně nainstalované verze firmwaru.
- Připojte USB flash disk do USB konektoru míchadla. Umístění USB konektorů najdete v oddíle "Přehled výrobku" na straně 44.
- 9. Zadejte přístupové heslo administrátora.



Obraz 114: Obrazovka Firmware

10. Stiskněte tlačítko Aktualizovat firmware.

UPOZORNĚNÍ Nedetekuje-li míchadlo žádný USB flash disk, zobrazí se výzva Připojte USB flash disk s firmwarem, který chcete nainstalovat, aby bylo možné spustit aktualizaci. Nyní připojte USB flash disk nebo zkontrolujte, zda je flash disk správně zastrčený do USB konektoru.

 Je-li k dispozici aktualizace firmwaru, zobrazí se na obrazovce Aktualizace firmwaru seznam znázorněný na Obraz 115. Stisknutím zvolte firmwarové položky, které chcete zaktualizovat.



Obraz 115: Výběr firmwarových položek pro aktualizaci

UPOZORNĚNÍ Firmwarové položky, jejichž verze jsou aktuální, jsou označené šedou barvou a není je možné zvolit. Jsou-li všechny verze firmwaru aktuální, zobrazí se hlášení V tuto chvíli nejsou k dispozici žádné nové aktualizace.

- 12. Stiskněte tlačítko Vybrat.
- Obsahuje-li flash disk několik verzí firmwaru, budete vyzváni k výběru jedné verze. Označte požadovanou verzi firmwaru a pak stiskněte tlačítko Vybrat.



Obraz 116: Výběr verze firmwaru pro instalaci

UPOZORNĚNÍ Nejste-li si jisti, kterou verzi máte nainstalovat, kontaktujte autorizovaného technika, od něhož jste obdrželi odkaz pro stažení.

 Zobrazí se seznam položek, které se mají zaktualizovat. Stiskněte tlačítko OK, aktualizace se zahájí.



Obraz 117: Seznam položek, které se mají zaktualizovat

UPOZORNĚNÍ Uživatelské rozhraní automaticky spravuje vzájemně závislé položky: Po označení firmwarové položky pro aktualizaci software automaticky přidá všechny další položky potřebné pro instalaci.

UPOZORNĚNÍ Označíte-li v seznamu na Obraz 115 volbu H.M.I., aby se provedla pouze aktualizace uživatelského rozhraní, nebude obrazovka uživatelského rozhraní po spuštění aktualizace po dobu až jedné minuty reagovat na další stisknutí aktivní oblasti ovládacího panelu. Míchadlo během této doby <u>nevypínejte</u>, aktualizace se jinak nedokončí. Po krátkém čase míchadlo znovu naběhne a je opět připravené k použití.

UPOZORNĚNÍ Označíte-li v seznamu na Obraz 115 dvě nebo tři volby, aby se provedla aktualizace <u>několika</u> firmwarových komponent současně, nesmí dojít k vypnutí míchadla během celého aktualizačního procesu.

15. Zobrazí se různá hlášení, která informují o průběhu instalace, jak je znázorněno na Obraz 118.



Obraz 118: Hlášení během aktualizace firmwaru

 Po zobrazení výzvy k restartování míchadla zařízení vypněte a znovu zapněte síťovým vypínačem.

4.8. Výměna základny

Základní pokyny pro montáž základny najdete v oddíle "3. 3. 1. Montáž základny" na straně 89. Kompletní seznam náhradních základen pro každé míchadlo najdete v oddíle "1. 2. 1. Základny" na straně 18.

UPOZORNĚNÍ Míchadla s regulací teploty je dovoleno provozovat jen se základnami předinstalovanými z výroby.

4.9. Servis

Společnost Thermo Fisher Scientific doporučuje nechat jednou ročně provést údržbu míchadla a příslušenství autorizovaným servisním technikem. Servisní technik kontroluje:

- elektrické instalace
- způsobilost místa instalace
- bezpečnostní systém
- použité příslušenství
- upevnění držáků baněk, základen a ostatního příslušenství k míchadlu

Před zahájením servisu je míchadlo a příslušenství nutné důkladně vyčistit a dekontaminovat, aby bylo možné provést úplnou a bezpečnou inspekci.

Pro tyto služby nabízí společnost Thermo Fisher Scientific možnost uzavření inspekčních a servisních smluv. Případné opravy jsou v rámci záručních podmínek prováděny zdarma, mimo rámec za úplatu. To platí jen v případě, že zásahy na míchadle byly prováděny výhradně servisními techniky společnosti Thermo Fisher Scientific.

4. 10. Odeslání a likvidace

VÝSTRAHA Poškození zdraví infekčními látkami. Při vyřazování míchadla a příslušenství z provozu pro následnou likvidaci je nutné celý systém vyčistit a v případě potřeby dezinfikovat nebo dekontaminovat. V případě pochybností kontaktujte zákaznický servis společnosti Thermo Fisher Scientific.

Při likvidaci míchadla dodržujte předpisy platné ve vaší zemi. Při likvidaci míchadla se obraťte na zákaznický servis společnosti Thermo Fisher Scientific. Kontaktní informace najdete na zadní straně tohoto návodu nebo na internetové stránce <u>www.thermofisher.com</u>.

Dodržte informace k přepravě a odeslání ("Přeprava" na straně 52 a "Odeslání" na straně 59).

5. Příčiny a odstraňování poruch

Č.	Označení	Řešení	Symbol
Chyba			
11999	Došlo k interní chybě	Došlo k interní chybě. Zařízení vypněte a znovu zapněte síťovým vypínačem. Zobrazuje-li se chybové hlášení i nadále, kontaktujte servisního technika.	
Výstražná	hlášení		
3	Vypršel interval pro kalibraci teploty. Uplynul 1 rok.	Vypršela roční platnost. Proveďte novou kalibraci teploty podle popisu v oddíle "4. 6. Kalibrace teploty" na straně 126.	CAL
4	Napájení obnoveno, aktivován automatický restart.	Při posledním zpracování došlo k výpadku elektrického proudu. Po obnově napájení se zpracování automaticky znovu zahájilo.	Ę
6	Výpadek senzorů teploty okolí.	Teplota naměřená senzorem teploty okolí je neplatná.	III
10	Otáčky ventilátoru 1 leží nad mezní hodnotou.	Otáčky ventilátoru 1 leží nad platnou mezí.	
11	Otáčky ventilátoru 2 leží nad mezní hodnotou.	Otáčky ventilátoru 2 leží nad platnou mezí.	•2
12	Otáčky ventilátoru 3 leží nad mezní hodnotou.	Otáčky ventilátoru 3 leží nad platnou mezí.	•2
13	Otáčky ventilátoru 4 leží nad mezní hodnotou.	Otáčky ventilátoru 4 leží nad platnou mezí.	•2
20	Kryt je již delší dobu otevřený.	Pro udržení teploty komory zavřete kryt míchadla.	4
Alarmová I	llášení		
1	Alarm nadměrné teploty.	Teplota komory překročila horní alarmovou hodnotu. Zkontrolujte vzorky, podmínky okolí a/nebo upravte nastavení.	
2	Alarm příliš nízké teploty.	Teplota komory překročila spodní alarmovou hodnotu. Zkontrolujte vzorky, podmínky okolí a/nebo upravte nastavení.	
3	Automatický restart po výpadku proudu se nezdařil.	Při posledním zpracování došlo k výpadku elektrického proudu. Nebylo možné úspěšně provést automatický restart.	Ċ
5	Příliš pomalá akcelerace pohonu. Není možné dosáhnout požadovaných otáček.	Nebylo možné dosáhnout požadovaných otáček. Zkontrolujte nastavení a/nebo náklad (příslušenství a vzorky) na základně.	\odot
7	Měření otáček ventilátoru 1 detekovalo neočekávaně klidový stav.	Zařízení vypněte a znovu zapněte. Zobrazuje-li se chybové hlášení i nadále, kontaktujte servisního technika.	•2
8	Měření otáček ventilátoru 2 detekovalo neočekávaně klidový stav.	Zařízení vypněte a znovu zapněte. Zobrazuje-li se chybové hlášení i nadále, kontaktujte servisního technika.	•2

Odeslání a likvidace

Č.	Označení	Řešení	Symbol
9	Měření otáček ventilátoru 3 detekovalo neočekávaně klidový stav.	Zařízení vypněte a znovu zapněte. Zobrazuje-li se chybové hlášení i nadále, kontaktujte servisního technika.	•2
10	Měření otáček ventilátoru 4 detekovalo neočekávaně klidový stav.	Zařízení vypněte a znovu zapněte. Zobrazuje-li se chybové hlášení i nadále, kontaktujte servisního technika.	•2
12	Chyba při rozběhu pohonu - chybí signál otáček.	Základna míchadla je zablokovaná. Zkontrolujte, že je okolo základny k dispozici dostatek místa, a/nebo zredukujte náklad (příslušenství a vzorky) na základně. Poté stiskněte na míchadle znovu tlačítko START. Zobrazuje-li se chybové hlášení i nadále, kontaktujte servisního technika.	\odot
24	Chyba měření otáček při běhu zařízení.	Detekovány nezvyklé otáčky. Zkontrolujte náklad (příslušenství a vzorky) a/nebo držáky baněk na základně. Poté stiskněte na míchadle znovu tlačítko START. Zobrazuje-li se chybové hlášení i nadále, kontaktujte servisního technika.	-
26	Měření otáček detekovalo při běhu neočekávaně klidový stav.	Zkontrolujte, že je okolo základny k dispozici dostatek místa, a/nebo snižte náklad (příslušenství a vzorky) na základně. Zkontrolujte, že nezareagovala pojistka na zadní straně míchadla ("Pojistka" na straně 47). Poté stiskněte na míchadle znovu tlačítko START. Zobrazuje-li se chybové hlášení i nadále, kontaktujte servisního technika.	\odot
82	Měření proudu motoru detekovalo přetižení.	Detekován nadměrný proud motoru. Nenakládejte/nevykládejte základnu při běhu zařízení. Snižte otáčky nebo upravte náklad (příslušenství a vzorky) na základně.	0
83	Měřené hodnoty proudu motoru leží mimo přípustný rozsah.	Detekován nadměrný proud motoru. Nenakládejte/nevykládejte základnu při běhu zařízení. Snižte otáčky nebo upravte náklad (příslušenství a vzorky) na základně.	— 0

Tabulka 59: Seznam chybových, výstražných a alarmových hlášení

UPOZORNĚNÍ Zobrazí-li se chybové hlášení, které není uvedené v této tabulce, kontaktujte servisního technika.

GPL (General Public License, obecná veřejná licence)

Část softwaru zařízení využívá otevřený software šířený pod GPL, LGPL nebo jinou open source licencí. Konkrétně se jedná o knihovny uvedené v tabulce níže. Zdrojový kód použitých knihoven (knihovny třetích stran) je možné získat u společnosti Thermo Fisher Scientific, pokud to příslušná licence dovoluje. Příslušné licenční podmínky použitého otevřeného softwaru jsou součástí poskytnutého balíku se zdrojovým kódem.

Knihovny třetích stran

Knihovna	Verze	Forma distribuce	Licence
Qt	5.8	BSP	LGPLv3
Log4Cplus	1.2.0	aplikace	veřejná licence Apache v2 / 2bodová BSD licence
boost	1.72.0	aplikace	licence Boost 1.0
json11	1.0.0	aplikace	licence MIT

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

Rejstřík

Α

Alarmová 69 Autoklávování 125 Automatický restart 73

B

Bezpečné otáčky 117

С

Čas 75 Časový režim 73 Čištění 123

D

Dekontaminace 124 Dezinfekce 124 Displej 73 Doba ztlumení 70 Dvoupatrové základny 91

Е

Ethernet 47

G

GPL (General Public License, obecná veřejná licence) 137 Grafické uživatelské rozhraní 60 Grafy 85

Н

Hlasitost 69

Instalace čtvercového držáku baněk 94 Instalace držáku pro dělicí nálevku 102 Instalace držáku pro mikrotitrační/DeepWell destičky 98 Instalace držáku stojanu na zkumavky s nastavitelným úhlem 97 Instalace nastavitelného držáku nádob 99 Instalace rozdělovače plynu 113 Instalace stojanu na zkumavky 96 Instalace stojanu pro kádinky 101 Instalace univerzálního odkládacího tácu 105

J

Jas 73 Jazyk 74

K

Kalibrace teploty 126 Klidový režim 75 Kontrola příslušenství 123 Kovové díly 123

L

Likvidace 134

М

Maximální otáčky 117 Místo instalace 50 Montáž základny 89

Ν

Nakládání 115 Nasazení držáků baněk a nádob 92 Nastavení 68 Název zařízení 76 Normy 43

0

Obsah dodávky 49 Obsluha 60 Odeslání 59, 134 Ρ

Péče 121 Plastové díly 123 Pojistka 47 Položení a použití adhezní podložky 106 Použití v souladu s určením 5, 115, 117 Požadovaná teplota 73 Požadované otáčky 73 Programy 78 Protokol událostí 84 Provozní parametry 73 Přehled výrobku 44 Přeprava 52 Příciny a odstraňování poruch 135 Přípojky 46 Příslušenství 17, 88

R

Region 76

S

Signální slova a symboly 6 Síťový konektor 47 Směrnice 43 Soubory a informace 77 Symboly použité v návodu 7

Т

Technické specifikace 11 Technické údaje 11

U

Údržba 121 Univerzální základny 90 USB 48

۷

Vkládání 116 Výstražná hlášení 69

Z

Základní nastavení 54 Základny 18

thermo scientific

Vyrobeno pro





Thermo Electron LED GmbH Zweigniederlassung Osterode Am Kalkberg, 37520 Osterode am Harz Germany Země původu: 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 je originální návod k použití. Tento návod k použití je překladem originálního návodu k použití.

thermofisher.com/shaker

© 2019–2023 Thermo Fisher Scientific Inc. Všechna práva vyhrazena.

Není-li výslovně uvedeno jinak, jsou všechny ochranné známky vlastnictvím společnosti Thermo Fisher Scientific Inc. a přidružených společností. Ne všechny výrobky jsou dostupné ve všech zemích. Přesnější informace získáte na požádání u místního prodejního partnera.

Obrázky publikované v tomto návodu jsou jen referenční. Zobrazená nastavení a jazyky se mohou lišit.

Austrálie +61 39757 4300

Rakousko +43 1 801 40 0

Belgie +32 53 73 42 41

Čína +800 810 5118 nebo +400 650 5118

Francie +33 2 2803 2180

Německo, bezplatná vnitrostátní linka 0800 1 536 376

Německo, mezinárodní linka +49 6184 90 6000 Indie +91 22 6716 2200

Itálie +39 02 95059 552

Japonsko +81 3 5826 1616

Nizozemí +31 76 579 55 55

Nový Zéland +64 9 980 6700

Severské země/Pobaltí/SNS země +358 10 329 2200

Rusko +7 812 703 42 15 Španělsko/Portugalsko +34 93 223 09 18

Švýcarsko +41 44 454 12 12

UK/Irsko +44 870 609 9203

USA/Kanada +1 866 984 3766

Ostatní asijské země +852 2885 4613

Ostatní země +49 6184 90 6000

