

Thermo Scientific Agitatori della serie Solaris

Istruzioni

70900193-е • 2024-11



Contenuto

Prefazione	9	. 5
Informazi	oni riguardanti questo manuale	. 5
Uso previ	sto	. 5
	Agitatore aperto	. 5
	Agitatore-incubatore (riscaldato) e agitatore refrigerato	_
	(raffreddato e riscaldato)	. 5
	Competenze richieste	. 5
Parole e	simboli di segnalazione	. 6
	Simboli utilizzati sull'agitatore e sugli accessori	. 6
	Simboli utilizzati nelle istruzioni	. 7
Avvertenz	ze di sicurezza	. 7
1. Specific	he tecniche	12
1. 1. Dat	i tecnici	12
	Agitatore Thermo Scientific Solaris 2000	.12
	Agitatore Thermo Scientific Solaris 4000	.13
	Agitatore Thermo Scientific Solaris 2000 I	.14
	Agitatore Thermo Scientific Solaris 4000 I	.15
	Agitatore Thermo Scientific Solaris 2000 R	.16
	Agitatore Thermo Scientific Solaris 4000 R	.17
1.2. Acc	essori	18
1. 2. 1.	Piattaforme	19
1. 2. 2.	Supporti porta matracci	20
1. 2. 3.	Supporti per micropiastra / piastra DeepWell	24
1. 2. 4.	Rack portaprovette	25
1. 2. 5.	Supporti ad angolo regolabile per rack portaprovette	29
1. 2. 6.	Supporti regolabili per recipienti	31
1. 2. 7.	Supporti per recipienti quadrati	33
1. 2. 8.	Rack porta becher Nalgene	35
1. 2. 9.	Supporto per imbuto separatore	39
1. 2. 10	. Supporto multiuso	40
1. 2. 11	. Tappetini antiscivolo	41
1. 2. 12	Accessori generali	43

1. 3.	Direttive e norme	14
1.4.	Informazioni generali sul prodotto	15
1.4	. 1. Solaris 2000	45
1.4	. 2. Solaris 4000	45
1.4	. 3. Solaris 2000 I / 2000 R	46
1.4	4. 4. Solaris 4000 I / 4000 R 4	46
1.4	. 5. Raccordi	47
2. Tras	porto e installazione	50
2.1.	Apertura dell'imballo	50
	Dotazione di fornitura	50
2. 2.	Luogo di installazione	51
2. 3.	Trasporto	53
	Manipolazione dell'agitatore.	53
2. 4.		54
2. 5.	Collegamento alla rete	54
2.6	Configurazione iniziale	55
2.0.		20
2. 7.		
2. 8.	Spedizione	50
3. Funz	zionamento6	61
3.1.	Accensione/spegnimento	61
3. 2.	Interfaccia utente grafica	51
3. 2	2. 1. Impostazione dei parametri operativi base per l'agitatore	62
3. 2	2. 2. Stato	67
3. 2	2. 3. Impostazioni	69
3. 2	2.4. Programmi	79
3. 3.	Accessori 8	39
3. 3	3. 1. Installazione della piattaforma	90
3. 3	8.2. Installazione dei supporti porta matracci e dei recipienti 9	93
3. 3	8.3. Installazione dei supporti quadrati	95
3. 3	8.4. Installazione dei rack portaprovette	97
3. 3	5. 5. Installazione dei supporti ad angolo regolabile per rack	20
port	taprovette	98

3. 3. 6. Installazione dei supporti per micropiastra / piastra	
	100
3. 3. 7. Installazione del supporti regolabili per recipienti	100
3. 3. 9. Installazione del supporto per imbuto separatore	102
3. 3. 10. Installazione del supporto multiuso	106
3. 3. 11. Installazione e uso del tappetino antiscivolo	107
3. 3. 12. Installazione del collettore del gas	114
3. 4. Caricamento e uso previsto	116
Caricamento	117
Uso previsto	118
4. Manutenzione e cura	122
4. 1. Principi fondamentali	123
Ispezione degli accessori	124
4. 2. Pulizia	124
Touch-screen	124
4. 3. Disinfezione	125
4. 4. Decontaminazione	125
4. 5. Autoclavaggio	126
4. 6. Calibrazione di temperatura	127
4. 6. 1. Buone pratiche per la calibrazione della temperatura .	127
4. 6. 2. Effettuazione della calibrazione di temperatura	127
4. 6. 3. Individuazione della temperatura calibrata	131
4. 7. Installazione del firmware	131
4. 8. Sostituzione della piattaforma	135
4.9. Manutenzione	135
4. 10. Spedizione e smaltimento	135
5. Eliminazione guasti	136
GPL (Licenza Pubblica Generica)	138
Indice analitico	139

Prefazione

Informazioni riguardanti questo manuale

Questo manuale di istruzioni descrive gli agitatori elencati di seguito e gli accessori compatibili («1. 2. Accessori» a pagina 18):

Agitatore	No. art.
Agitatore Thermo Scientific Solaris 2000, 100–240 V ± 10 %, 50 / 60 Hz	SK2000
Agitatore Thermo Scientific Solaris 4000, 100–240 V ± 10 %, 50 / 60 Hz	SK4000
Agitatore Thermo Scientific Solaris 2000 I, 100–120, 200–240 V ± 10 %, 50 / 60 Hz	SK2001
Agitatore Thermo Scientific Solaris 4000 I, 100–120, 200–240 V ± 10 %, 50 / 60 Hz	SK4001
Agitatore Thermo Scientific Solaris 2000 R, 100–240 V ± 10 %, 50 / 60 Hz	SK2002
Agitatore Thermo Scientific Solaris 4000 R, 100–240 V ±10 %, 50 / 60 Hz	SK4002

Anziché indicare il nome completo del prodotto, le descrizioni e istruzioni contenute in questo manuale fanno riferimento semplicemente alle singole classi di agitatori:

- Agitatore aperto
- Agitatore-incubatore (riscaldato)
- Agitatore refrigerato (raffreddato e riscaldato)

Uso previsto

Agitatore aperto

Questo dispositivo viene utilizzato nel campo della ricerca e della produzione ed è destinato all'agitazione di soluzioni in un'orbita 2D, controllando al contempo la velocità e il tempo nelle applicazioni.

Agitatore-incubatore (riscaldato) e agitatore refrigerato

(raffreddato e riscaldato)

Questo dispositivo viene utilizzato nel campo della ricerca e della produzione ed è destinato all'agitazione di soluzioni in un'orbita 2D, controllando al contempo la velocità, la durata e la temperatura nelle applicazioni.

Competenze richieste

Questo agitatore deve essere utilizzato solo da personale istruito.

Parole e simboli di segnalazione

Parola di segnalazione	Livello di pericolo
AVVERTENZA	Sta ad indicare una situazione di pericolo che causerà la morte o lesioni gravi se non viene evitata.
ATTENZIONE	Sta ad indicare una situazione di pericolo che causerà lesioni minori o moderate, se non viene evitata.
AVVISO	Sta ad indicare informazioni considerate importanti, ma non legate a pericoli.

Simboli utilizzati sull'agitatore e sugli accessori

Rispettare le indicazioni nel manuale per non mettere in pericolo se stessi e l'ambiente.

(Leggere le istruzioni per l'uso
	Staccare la spina
	Pericolo generale
	Pericolo causato da superficie molto calda
	Rischio di lesioni alle dita o alle mani causate da parti meccaniche che si chiudono all'improvviso

Simboli utilizzati nelle istruzioni

Rispettare le indicazioni nel manuale per non mettere in pericolo se stessi e l'ambiente.

Pericolo generale	<u>A</u>	Pericolo elettrico
Rischio biologico		Rischio di taglio
Pericolo causato da materiali infiammabili	i	Sta ad indicare informazioni considerate importanti, ma non legate a pericoli.
Rischio di lesioni alle dita o alle mani causate da parti meccaniche che si chiudono all'improvviso.		

Avvertenze di sicurezza



La mancata osservanza di queste avvertenze di sicurezza può portare a situazioni pericolose che, se non evitate, possono provocare la morte o gravi lesioni.

AVVERTENZA

- Rispettare le informazioni per la sicurezza.
- Utilizzare l'agitatore solo in conformità all'uso previsto. Un uso improprio può causare danni, contaminazioni e lesioni anche mortali.
- L'agitatore deve essere utilizzato solo da personale adeguatamente istruito.
- Il gestore è obbligato a garantire che vengano indossati i dispositivi di protezione individuale idonei. L'operatore dovrebbe conoscere il manuale di biosicurezza nei laboratori «Laboratory Biosafety Manual» (dell'Organizzazione Mondiale della Sanità) riconosciuto a livello internazionale, come anche le raccomandazioni nazionali vigenti in materia.



Danni causati da un'alimentazione elettrica inadatta.

Assicurarsi che l'agitatore sia collegato sempre solo a prese correttamente messe a terra.



Rischi provocati dalla manipolazione di sostanze pericolose.

AVVERTENZA

Quando si lavora con campioni corrosivi (soluzioni saline, acidi, basi) gli accessori e l'agitatore devono essere puliti accuratamente.

- L'agitatore non è né inerte né antideflagrante. Non utilizzare mai l'agitatore in un ambiente a rischio di deflagrazione.
- Non agitare materiali tossici o radioattivi o microorganismi patogeni senza misure di sicurezza idonee.
- Per l'agitazione di materiali pericolosi attenersi al manuale di biosicurezza nei laboratori («Laboratory Biosafety Manual») dell'Organizzazione Mondiale della Sanità (OMS) e alle norme nazionali vigenti in materia. Se vengono agitate prove microbiologiche del gruppo di rischio II (secondo il manuale «Laboratory Biosafety Manual» dell'organizzazione di salute mondiale WHO) devono essere impiegate chiusure bio-aerosol con tenuta aerosol. Sul sito dell'Organizzazione Mondiale della Sanità OMS (www.who.int); (WHO, World Health Organisation) cercare il manuale «Laboratory Biosafety Manual». Con materiali con un gruppo di rischio più alto deve essere previsto più di un provvedimento di protezione.
- Se l'agitatore o i suoi componenti sono stati contaminati da sostanze tossiche o patogene, è necessario adottare misure di disinfezione e decontaminazione idonee («Decontaminazione» a pagina 125; «Disinfezione» a pagina 125).
- Al verificarsi di una situazione di pericolo, disinserire l'alimentazione dell'agitatore e abbandonare subito l'ambiente circostante l'agitatore.



Danni alla salute provocati da sostanze infettive.

AVVERTENZA

Qualora liquidi o altri materiali dovessero essere versati accidentalmente sotto la piattaforma, spegnere subito l'agitatore, staccare la spina e rimuovere la piattaforma («Piattaforme» a pagina 19).

Eliminare il materiale versato seguendo le regolari procedure adottate nel laboratorio. Utilizzare i dispositivi di protezione individuale idonei.



Danni alla salute causati dall'agitazione di sostanze o materiali esplosivi o infiammabili. Non agitare sostanze o materiali esplosivi o infiammabili.

AVVERTENZA



ATTENZIONE

Ustioni causate da superfici calde.

In un agitatore a termoregolazione le superfici, come ad esempio la piattaforma e gli accessori, possono riscaldarsi fortemente sotto la cappa. Non toccare le superfici calde. Aspettare che le superfici calde si siano raffreddate.



Ferite da taglio causate da frammenti di vetro.

ATTENZIONE

Una piattaforma o accessori, come ad es. supporti porta beute, allentati possono comportare frammenti di vetro a seguito di recipienti che cadono dall'agitatore. Assicurarsi che la piattaforma e gli accessori siano montati accuratamente, con l'ausilio degli attrezzi e delle viti corretti. Assicurarsi che i supporti porta beute siano in grado di sostenere il carico del recipiente e del campione alla velocità scelta. I supporti porta beute con carichi configurati individualmente potrebbero avere una velocità massima inferiore ai limiti di stabilità dell'agitatore. Consultare «3. 4. Caricamento e uso previsto» a pagina 116.

Badare a rumori insoliti. Questi potrebbero indicare una piattaforma o accessori allentati.



Danni all'agitatore o malfunzionamento causati da un touch-screen danneggiato.

ATTENZIONE

- Non usare l'agitatore con il touch-screen danneggiato. Spegnere l'agitatore. Staccare la spina. Fare sostituire
- il touch-screen da un tecnico autorizzato del servizio di assistenza.



Rischio biologico dovuto a recipienti rotti o che perdono.

Accessori non installati correttamente possono comportare la fuoriuscita dei campioni.

ATTENZIONE

- Assicurarsi che gli accessori siano installati accuratamente con l'ausilio degli attrezzi e delle viti corretti.
- Assicurarsi che gli accessori siano adatti alla piattaforma.
- Utilizzare i recipienti sempre con gli accessori adatti alla rispettiva dimensione.
- I recipienti devono essere intatti e installati accuratamente.



La sicurezza può essere pregiudicata da un caricamento non corretto e da accessori difettosi.

ATTENZIONE

- Assicurarsi sempre che il carico (accessori e campioni) sia distribuito il più uniformemente possibile, specialmente quando si lavora con una piattaforma a due livelli.
- Non usare accessori che presentano tracce di corrosione o incrinature. Per maggiori informazioni contattare il servizio di assistenza ai clienti.
- Utilizzare solo agitatori che sono stati caricati correttamente.
- Non sovraccaricare mai l'agitatore.
- Assicurarsi che gli accessori siano stati installati correttamente prima di usare l'agitatore. Attenersi alle istruzioni riportate al capitolo «Accessori» a pagina 89.



Lesioni causate dalla mancata osservanza dei principi fondamentali.

ATTENZIONE

- Non usare mai l'agitatore senza piattaforma installata correttamente.
- Non utilizzare mai l'agitatore se parti del corpo sono danneggiate o sono state rimosse.
- Non usare mai un agitatore con la molla della cappa difettosa. Una molla difettosa non è in grado di mantenere la cappa aperta in sicurezza nella posizione completamente aperta.
- Non spostare l'agitatore mentre è in funzione.
- Non appoggiarsi sull'agitatore.
- Non caricare o scaricare mai l'agitatore prima che abbia raggiunto definitivamente la condizione di fermo e questa sia confermata sul touch-screen.
- Non mettere mai oggetti sull'agitatore mentre questo è in funzione.
- Non toccare la piattaforma o gli accessori sull'agitatore mentre questo è in funzione.
- Il corpo dell'agitatore non deve essere aperto dall'operatore.



La protezione può essere pregiudicata da accessori incompatibili.

AVVISO

Utilizzare esclusivamente accessori omologati da Thermo Fisher Scientific per quest'agitatore. Per gli elenchi aggiornati visitare <u>www.thermofisher.com</u>.



AVVISO

Per arrestare l'agitatore:

Premere il tasto STOP. Spegnere l'agitatore azionando l'interruttore generale. Staccare la spina di alimentazione elettrica. In caso di emergenza disconnettere l'alimentazione elettrica.

1. Specifiche tecniche

1.1. Dati tecnici

Agitatore Thermo Scientific Solaris 2000					
Gamma di velocità	15–525 giri/min				
Tempo di funzionamento	99 ore 59 min (incremento di 1 min) o funzionamento continuo				
Livello sonoro alla velocità massima	50 dB (A) (a 1 m di dista	anza dal dispositivo e a un	'altezza di 1,6 m)		
Carico massimo (compr. piattaforma, accessori e campioni)	25 kg (55 lbs)				
Alimentazione elettrica	100–240 V ±10 %, 50 /	/ 60 Hz			
Potenza assorbita	60 W				
Condizioni ambientali					
Per lo stoccaggio e il trasporto	Temperatura: da -10 °C Umidità: da 15 % a 85 9	a 55 °C %			
Per l'esercizio	Utilizzo in ambiente chiusi Altitudini fino a 3 000 metri sopra il livello del mare Temperatura: da 5 °C a 40 °C Umidità relativa massima dell'80 % per temperature sino a 31 °C, decrescente linearmente sino ad umidità relativa 50 % a 40 °C				
Grado di inquinamento	2				
Categoria di sovratensione	Ш				
IP	20				
Interfacce	USB	Ethernet			
	2x USB-A 2.0	RJ45			
Dimensioni	Lunghezza	Larghezza	Altezza		
	47 cm (18,5 in)	37 cm (14,5 in)	15 cm (5,5 in)		
Peso	20.9 kg (46.0 lbs)				

Tabella 1: Solaris 2000 - Dati tecnici

Agitatore Thermo Scientific Solaris 4000

Gamma di velocità	15–525 giri/min				
Tempo di funzionamento	99 ore 59 min (incremento di 1 min) o funzionamento continuo				
Livello sonoro alla velocità massima	54 dB (A) (a 1 m di dista	54 dB (A) (a 1 m di distanza dal dispositivo e a un'altezza di 1,6 m)			
Carico massimo (compr. piattaforma, accessori e campioni)	43 kg (95 lbs)	43 kg (95 lbs)			
Alimentazione elettrica	100–240 V ±10 %, 50 /	60 Hz			
Potenza assorbita	80 W				
Condizioni ambientali					
Per lo stoccaggio e il trasporto	Temperatura: da -10 °C Umidità: da 15 % a 85 9	a 55 °C 6			
Per l'esercizio	Utilizzo in ambiente chiusi Altitudini fino a 3 000 metri sopra il livello del mare Temperatura: da 5 °C a 40 °C Umidità relativa massima dell'80 % per temperature sino a 31 °C, decrescente linearmente sino ad umidità relativa 50 % a 40 °C				
Grado di inquinamento	2				
Categoria di sovratensione	I				
IP	20				
Interfacce	USB	Ethernet			
	2x USB-A 2.0	RJ45			
Dimensioni	Lunghezza	Larghezza	Altezza		
	65 cm (25,6 in)	58 cm (22,8 in)	18 cm (7 in)		
Peso	75,1 kg (165,5 lbs)				

Tabella 2: Solaris 4000 - Dati tecnici

Agitatore Thermo Scientific Solaris 2000 I

Gamma di velocità	15–525 giri/min		
Range di impostazione temperatura	30-60 °C		
Stabilità della temperatura nel matraccio a 37 °C	Temperatura ambiente 23 °C; per 1 ora con temperatura stabile nella camera \pm 0,1 °C		
Uniformità della temperatura nel matraccio a 37 °C	Temperatura ambiente 23 ± 0,5 °C	°C; per 1 ora con temperat	ura stabile nella camera
Tempo di funzionamento	99 ore 59 min (incremen	to di 1 min) o funzionamen	to continuo
Livello sonoro alla velocità massima	52 dB (A) (a 1 m di dista	nza dal dispositivo e a un'a	litezza di 1,6 m)
Carico massimo (compr. piattaforma, accessori e campioni)	16 kg (35 lbs), compresa piattaforma 1	1 x 14 da 2,3 kg (5,1 lbs)	
Alimentazione elettrica	100–120, 200–240 V ±1	10 %, 50 / 60 Hz	
Potenza assorbita	900 W		
Condizioni ambientali			
Per lo stoccaggio e il trasporto	Temperatura: da -10 °C a 55 °C Umidità: da 15 % a 85 %		
Per l'esercizio	Utilizzo in ambiente chiusi Altitudini fino a 3 000 metri sopra il livello del mare Temperatura: da 5 °C a 40 °C Umidità relativa massima dell'80 % per temperature sino a 31 °C, decrescente linearmente sino ad umidità relativa 50 % a 40 °C		
Grado di inquinamento	2		
Categoria di sovratensione			
IP	20		
Interfacce	USB	Ethernet	
	2x USB-A 2.0	RJ45	
Dimensioni	Lunghezza	Larghezza	Altezza
	70 cm (27,6 in)	36 cm (14,2 in)	46 cm (18,1 in)
Con cappa aperta			79 cm (30,8 in)
Peso	44,9 kg (99 lbs)		

Tabella 3: Solaris 2000 I - Dati tecnici

Agitatore Thermo Scientific Solaris 4000 I

Gamma di velocità	15–525 giri/min		
Range di impostazione temperatura	30−60 °C		
Stabilità della temperatura nel matraccio a 37 °C	Temperatura ambiente 23 °C; per 1 ora con temperatura stabile nella camera \pm 0,1 °C		
Uniformità della temperatura nel matraccio a 37 °C	Temperatura ambiente 23 ± 0,5 °C	°C; per 1 ora con temperati	ura stabile nella camera
Tempo di funzionamento	99 ore 59 min (increment	to di 1 min) o funzionamen	to continuo
Livello sonoro alla velocità massima	52 dB (A) (a 1 m di distar	nza dal dispositivo e a un'a	iltezza di 1,6 m)
Carico massimo (compr. piattaforma, accessori e campioni)	32 kg (70,5 lbs), compresa piattaforma 18	3 x 18 da 4,3 kg (9,5 lbs)	
Alimentazione elettrica	100–120, 200–240 V ±1	0 %, 50 / 60 Hz	
Potenza assorbita	900 W		
Condizioni ambientali			
Per lo stoccaggio e il trasporto	Temperatura: da -10 °C a Umidità: da 15 % a 85 %	a 55 °C	
Per l'esercizio	Utilizzo in ambiente chiusi Altitudini fino a 3 000 metri sopra il livello del mare Temperatura: da 5 °C a 40 °C Umidità relativa massima dell'80 % per temperature sino a 31 °C, decrescente linearmente sino ad umidità relativa 50 % a 40 °C		
Grado di inquinamento	2		
Categoria di sovratensione	1		
IP	20		
Interfacce	USB	Ethernet	
	2x USB-A 2.0	RJ45	
Dimensioni	Lunghezza	Larghezza	Altezza
	77 cm (30,3 in)	57 cm (22,4 in)	55 cm (21,7 in)
Con cappa aperta			94 cm (36,7 in)

Tabella 4: Solaris 4000 I - Dati tecnici

Peso

68,1 kg (150,1 lbs)

Agitatore Thermo Scientific Solaris 2000 R

Gamma di velocità	15–525 giri/min		
Range di impostazione temperatura	5–60 °C		
Stabilità della temperatura nel matraccio a 37 °C	Temperatura ambiente 23 °C; per 1 ora con temperatura stabile nella camera \pm 0,1 °C		
Uniformità della temperatura nel matraccio a 37 °C	Temperatura ambiente 23 ± 0,5 °C	°C; per 1 ora con temperat	ura stabile nella camera
Tempo di funzionamento	99 ore 59 min (incremen	to di 1 min) o funzionamen	to continuo
Livello sonoro alla velocità massima	56 dB (A) (a 1 m di dista	nza dal dispositivo e a un'a	iltezza di 1,6 m)
Carico massimo (compr. piattaforma, accessori e campioni)	16 kg (35 lbs), compresa piattaforma 1	1 x 14 da 2,3 kg (5,1 lbs)	
Alimentazione elettrica	100–240 V ±10 %, 50 /	60 Hz	
Potenza assorbita	350 W		
Condizioni ambientali			
Per lo stoccaggio e il trasporto	Temperatura: da -10 °C a 55 °C Umidità: da 15 % a 85 %		
Per l'esercizio	Utilizzo in ambiente chiusi Altitudini fino a 3 000 metri sopra il livello del mare Temperatura: da 5 °C a 40 °C Umidità relativa massima dell'80 % per temperature sino a 31 °C, decrescente linearmente sino ad umidità relativa 50 % a 40 °C		
Grado di inquinamento	2		
Categoria di sovratensione	11		
IP	20		
Interfacce	USB	Ethernet	
	2x USB-A 2.0	RJ45	
Dimensioni	Lunghezza	Larghezza	Altezza
	70 cm (27,6 in)	36 cm (14,2 in)	46 cm (18,1 in)
Con cappa aperta			79 cm (30,8 in)
Peso	47,2 kg (104,1 lbs)		

Tabella 5: Solaris 2000 R - Dati tecnici

Agitatore Thermo Scientific Solaris 4000 R

Gamma di velocità	15–525 giri/min			
Range di impostazione temperatura	4–60 °C			
Stabilità della temperatura nel matraccio a 37 °C	Temperatura ambiente 23 ° ± 0,1 °C	°C; per 1 ora con temperatu	ura stabile nella camera	
Uniformità della temperatura nel matraccio a 37 °C	Temperatura ambiente 23 °C; per 1 ora con temperatura stabile nella camera \pm 0,5 °C			
Tempo di funzionamento	99 ore 59 min (increment	o di 1 min) o funzionament	to continuo	
Livello sonoro alla velocità massima	56 dB (A) (a 1 m di distan	za dal dispositivo e a un'a	ltezza di 1,6 m)	
Carico massimo (compr. piattaforma, accessori e campioni)	32 kg (70,5 lbs), compresa piattaforma 18	3 x 18 da 4,3 kg (9,5 lbs)		
Alimentazione elettrica	100–240 V ±10 %, 50 / 6	60 Hz		
Potenza assorbita	750 W			
Condizioni ambientali				
Per lo stoccaggio e il trasporto	Temperatura: da -10 °C a Umidità: da 15 % a 85 %	.55 ℃		
Per l'esercizio	Utilizzo in ambiente chiusi Attitudini fino a 3 000 metri sopra il livello del mare Temperatura: da 5 °C a 40 °C Umidità relativa massima dell'80 % per temperature sino a 31 °C, decrescente linearmente sino ad umidità relativa 50 % a 40 °C			
Grado di inquinamento	2			
Categoria di sovratensione	II			
P	20			
Interfacce	USB	Ethernet		
	2x USB-A 2.0	RJ45		
Dimensioni	Lunghezza	Larghezza	Altezza	
	77 cm (30,3 in)	57 cm (22,4 in)	55 cm (21,7 in)	
Con cappa aperta			94 cm (36,7 in)	
Peso	74,9 kg (165,1 lbs)			

Tabella 6: Solaris 4000 R - Dati tecnici

Accessori

1.2. Accessori



La protezione può essere pregiudicata da accessori incompatibili.

ATTENZIONE Utilizzare esclusivamente accessori omologati da Thermo Fisher Scientific per quest'agitatore.

Per gli elenchi aggiornati visitare <u>www.thermofisher.com</u>.

Accessori

1.2.1. Piattaforme

Piattaforma		
Solaris 2000	Peso	
Piattaforma universale Thermo Scientific Solaris 12x14	2,7 kg (5,9 lbs)	SK1214
Piattaforma universale a due livelli Thermo Scientific Solaris 12x14	6,4 kg (14,1 lbs)	SK1214D
Kit di upgrade Thermo Scientific Solaris 12x14 piattaforma universale a due livelli		SK1214DK
Piattaforma universale Thermo Scientific Solaris 18x18	4,3 kg (9,4 lbs)	SK1818
Plattaforma universale a due livelli Thermo Scientific Solaris 18x18	9,7 kg (21,3 lbs)	SK1818D
Kit di upgrade Thermo Scientific Solaris 18x18 piattaforma universale a due livelli		SK1818DK
Piattaforma universale Thermo Scientific Solaris 18x24	6,0 kg (13,2 lbs)	SK1824
Solaris 4000		
Plattaforma universale Thermo Scientific Solaris 18x30	7,5 kg (16,5 lbs)	SK1830
Piattaforma universale a due livelli Thermo Scientific Solaris 18x30	16,3 kg (35,9 lbs)	SK1830D
Kit di upgrade Thermo Scientific Solaris 18x30 piattaforma universale a due livelli		SK1830DK
Piattaforma universale Thermo Scientific Solaris 36x24	12,0 kg (26,4 lbs)	SK3624
Solaris 2000 I / 2000 R		
Piattaforma universale Thermo Scientific Solaris 11x14	2,3 kg (5,1 lbs)	SK1114
Solaris 4000 I / 4000 R		
Piattaforma universale Thermo Scientific Solaris 18x18	4,3 kg (9,5 lbs)	SK1818
Kit di ricambi e accessori		
Kit di ricambi per supporto (viti)		SK1001
Kit di ricambi per piattaforma, piccolo (viti per piattaforma, attrezzo)		SK0100
Kit di ricambi per piattaforma, grande (viti per piattaforma, attrezzo)		SK0101
Cacciavite per accessori		75004131

Tabella 7: Piattaforme disponibili

1. 2. 2. Supporti porta matracci



Per piattaforme Solaris 2000

	No. art.	Universale			A due livelli		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18	
Beuta da 10 ml	30150BI	72	113	157	140	226	
Beuta da 25 ml	30151	42	64	80	80	124	
Beuta da 50 ml	30152BI	42	64	80	80	124	
Beuta da 125 ml	30153	15	32	40	30	62	
Beuta da 250 ml	30154BI	9	16	24	16	32	
Beuta da 300 ml	30155	9	16	20	16	32	
Beuta da 500 ml	30156BI	9	16	20	16	32	
Beuta da 1 l	30157BI	4	9	11	8	16	
Beuta da 2 l	30158	3	5	6	-	-	
Beuta da 4 l	30159	1	4	4	-	-	
Beuta da 5 l	30159B	1	2	4	-	-	
Beuta da 6 l	30160	1	2	2	-	-	
Beuta di Fembach da 2800 ml	30162	1	4	4	-	-	
Beuta per colture da 2500 ml basso	30161	1	2	2	-	-	

 Tabella 8: Supporti porta matracci disponibili per piattaforme Solaris 2000

Per piattaforme Solaris 4000

	No. art.	Universale		A due livelli
		18 x 30	36 x 24	18 x 30
Beuta da 10 ml	30150BI	203	187	402
Beuta da 25 ml	30151	112	187	220
Beuta da 50 ml	30152BI	112	187	220
Beuta da 125 ml	30153	46	83	92
Beuta da 250 ml	30154BI	34	40	68
Beuta da 300 ml	30155	28	40	56
Beuta da 500 ml	30156BI	28	40	56
Beuta da 1 l	30157BI	14	20	28
Beuta da 2 l	30158	8	11	16
Beuta da 4 l	30159	6	8	-
Beuta da 5 l	30159B	5	8	-
Beuta da 6 I	30160	3	6	-
Beuta di Fembach da 2800 ml	30162	6	8	12
Beuta per colture da 2500 ml basso	30161	3	6	6

 Tabella 9:
 Supporti porta matracci disponibili per piattaforme Solaris 4000

Per piattaforme Solaris 2000 I / 2000 R

	No. art.	Universale
		11 x 14
Beuta da 10 ml	30150BI	59
Beuta da 25 ml	30151	35
Beuta da 50 ml	30152BI	35
Beuta da 125 ml	30153	15
Beuta da 250 ml	30154BI	7
Beuta da 300 ml	30155	7
Beuta da 500 ml	30156BI	7
Beuta da 1 l	30157BI	4
Beuta da 2 l	30158	2
Beuta da 4 l	30159	-
Beuta da 5 l	30159B	-
Beuta da 6 l	30160	-
Beuta di Fembach da 2800 ml	30162	1
Beuta per colture da 2500 ml basso	30161	1

Tabella 10: Supporti porta matracci disponibili per piattaforme Solaris 2000 I / 2000 R

Per piattaforme Solaris 4000 I / 4000 R

	No. art.	Universale
		18 x 18
Beuta da 10 ml	30150BI	113
Beuta da 25 ml	30151	64
Beuta da 50 ml	30152BI	64
Beuta da 125 ml	30153	32
Beuta da 250 ml	30154BI	16
Beuta da 300 ml	30155	16
Beuta da 500 ml	30156BI	16
Beuta da 1 l	30157BI	9
Beuta da 2 l	30158	6
Beuta da 4 l	30159	-
Beuta da 5 l	30159B	-
Beuta da 6 l	30160	-
Beuta di Fernbach da 2800 ml	30162	4
Beuta per colture da 2500 ml basso	30161	1

Tabella 11: Supporti porta matracci disponibili per piattaforme Solaris 4000 I / 4000 R

1. 2. 3. Supporti per micropiastra / piastra DeepWell



Per piattaforme Solaris 2000

	No. art.	Universale		A due livelli		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Micropiastra / piastra Deepwell	30175	7	12	16	14	24

Tabella 12: Capacità dei supporti per micropiastra / piastra DeepWell per piattaforme Solaris 2000

Per piattaforme Solaris 2000 I / 2000 R

	No. art.	Universale
Micropiastra / piastra Deepwell	30175	6

Tabella 13: Capacità dei supporti per micropiastra / piastra DeepWell per piattaforme Solaris2000 I / 2000 R

Per piattaforme Solaris 4000

	No. art.	Universale		A due livelli	
		18 x 30	36 x 24	18 x 30	
Micropiastra / piastra Deepwell	30175	23	36	42	

Tabella 14: Capacità dei supporti per micropiastra / piastra DeepWell per piattaforme Solaris 4000

Per piattaforme Solaris 4000 I / 4000 R

	No. art.	Universale
		18 x 18
Micropiastra / piastra Deepwell	30175	10

Tabella 15: Capacità dei supporti per micropiastra / piastra DeepWell per piattaforme Solaris4000 I / 4000 R

1. 2. 4. Rack portaprovette

Per piattaforme Solaris 2000



	No. art.	Universale		e	A due livelli	
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Formato ridotto						
10–13 mm, rosso, disposizione 6 x 6	30181	8	12	15	14	22
14–16 mm, arancione, disposizione 6 x 6	30183	5	9	11	9	16
17–20 mm, bianco, disposizione 4 x 5	30185	7	11	14	12	20
21–25 mm, azzurro, disposizione 4 x 4	30187	6	9	11	10	16
26–30 mm, verde, disposizione 3 x 3	30189	6	9	12	11	17
Microcentrifuga, 1,5 ml, azzurro, disposizione 4 x 6	30191	6	10	13	11	19
Formato normale						
10–13 mm, disposizione 6 x 12	30180BI	3	7	9	6	12
14–16 mm, disposizione 6 x 12	30182	3	4	6	6	8
17–20 mm, disposizione 4 x 10	30184	3	5	7	6	8
21–25 mm, disposizione 4 x 10	30186	2	3	5	4	6
26–30 mm, disposizione 3 x 8	30188	3	4	5	5	8
Microcentrifuga, 1,5 ml, disposizione 8 x 12	30190	З	4	7	6	8

Tabella 16: Rack portaprovette disponibili per piattaforme Solaris 2000

Accessori

Per piattaforme Solaris 4000

	No. art.	Unive	Universale	
		18 x 30	36 x 24	18 x 30
Formato ridotto				
10–13 mm, rosso, disposizione 6 x 6	30181	21	32	40
14–16 mm, arancione, disposizione 6 x 6	30183	13	20	24
17–20 mm, bianco, disposizione 4 x 5	30185	18	20	34
21–25 mm, azzurro, disposizione 4 x 4	30187	14	22	26
26–30 mm, verde, disposizione 3 x 3	30189	15	24	30
Microcentrifuga, 1,5 ml, azzurro, disposizione 4 x 6	30191	18	24	34
Formato normale				
10–13 mm, disposizione 6 x 12	30180BI	13	20	18
14–16 mm, disposizione 6 x 12	30182	9	12	16
17–20 mm, disposizione 4 x 10	30184	9	15	18
21–25 mm, disposizione 4 x 10	30186	6	9	12
26–30 mm, disposizione 3 x 8	30188	7	10	13
Microcentrifuga, 1,5 ml, disposizione 8 x 12	30190	9	12	17

Tabella 17: Rack portaprovette disponibili per piattaforme Solaris 4000

Per piattaforme Solaris 2000 I / 2000 R

	No. art.	Universale
Formato ridotto		
10–13 mm, rosso, disposizione 6 x 6	30181	6
14–16 mm, arancione, disposizione 6 x 6	30183	2
17–20 mm, bianco, disposizione 4 x 5	30185	4
21–25 mm, azzurro, disposizione 4 x 4	30187	2
26–30 mm, verde, disposizione 3 x 3	30189	4
Microcentrifuga, 1,5 ml, azzurro, disposizione 4 x 6	30191	3
Formato normale		
10–13 mm, disposizione 6 x 12	30180BI	3
14–16 mm, disposizione 6 x 12	30182	2
17–20 mm, disposizione 4 x 10	30184	2
21–25 mm, disposizione 4 x 10	30186	1
26–30 mm, disposizione 3 x 8	30188	2
Microcentrifuga, 1,5 ml, disposizione 8 x 12	30190	2

Tabella 18: Rack portaprovette disponibili per piattaforme Solaris 2000 I / 2000 R

Per piattaforme Solaris 4000 I / 4000 R

	No. art.	Universale
		18 x 18
Formato ridotto		
10–13 mm, rosso, disposizione 6 x 6	30181	12
14–16 mm, arancione, disposizione 6 x 6	30183	8
17–20 mm, bianco, disposizione 4 x 5	30185	10
21–25 mm, azzurro, disposizione 4 x 4	30187	8
26–30 mm, verde, disposizione 3 x 3	30189	9
Microcentrifuga, 1,5 ml, azzurro, disposizione 4 x 6	30191	8
Formato normale		
10–13 mm, disposizione 6 x 12	30180BI	7
14–16 mm, disposizione 6 x 12	30182	4
17–20 mm, disposizione 4 x 10	30184	5
21–25 mm, disposizione 4 x 10	30186	3
26–30 mm, disposizione 3 x 8	30188	4
Microcentrifuga, 1,5 ml, disposizione 8 x 12	30190	5

Tabella 19: Rack portaprovette disponibili per piattaforme Solaris 4000 I / 4000 R

1. 2. 5. Supporti ad angolo regolabile per rack portaprovette



Per piattaforme Solaris 2000

	No. art.	Universale			A due livelli		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18	
10-13 mm, 72 posti	236090	3	4	6	6	8	
16-20 mm, 40 posti	236091	3	4	6	5	8	
21-25 mm, 40 posti	236092	3	3	4	5	6	
26-30 mm, 24 posti	236093	3	З	4	5	6	

 Tabella 20:
 Supporti ad angolo regolabile per rack portaprovette disponibili per piattaforme

 Solaris 2000
 Solaris 2000

Per piattaforme Solaris 4000

	No. art.	Univ	ersale	A due livelli		
		18 x 30	36 x 24	18 x 30		
10-13 mm, 72 posti	236090	7	14	14		
16-20 mm, 40 posti	236091	7	14	14		
21-25 mm, 40 posti	236092	5	9	10		
26-30 mm, 24 posti	236093	6	9	12		

 Tabella 21: Supporti ad angolo regolabile per rack portaprovette disponibili per piattaforme
 Solaris 4000

Per piattaforme Solaris 2000 I / 2000 R

	No. art.	Universale
		11 x 14
10-13 mm, 72 posti	236090	2
16-20 mm, 40 posti	236091	2
21-25 mm, 40 posti	236092	1
26-30 mm, 24 posti	236093	1

Tabella 22: Supporti ad angolo regolabile per rack portaprovette disponibili per piattaformeSolaris 2000 I / 2000 R

Per piattaforme Solaris 4000 I / 4000 R

	No. art.	Universale
		18 x 18
10-13 mm, 72 posti	236090	4
16-20 mm, 40 posti	236091	3
21-25 mm, 40 posti	236092	3
26-30 mm, 24 posti	236093	3

Tabella 23: Supporti ad angolo regolabile per rack portaprovette disponibili per piattaforme Solaris 4000 I / 4000 R

1. 2. 6. Supporti regolabili per recipienti

Per piattaforme Solaris 2000



	No. art.	Universale			A due livelli		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18	
Supporto per recipiente singolo regolabile in altezza 11"x14"	75004104	-	-	-	-	-	
Supporto per recipiente a doppia regolazione in altezza 12"x14"	75004102	1	1	1	-	2	
Supporto per recipiente a doppia regolazione in altezza 18"x18"	75004103	-	1	1	-	-	
Supporto regolabile per recipiente singolo	75004101	4	4	6	6	8	

Tabella 24: Supporti regolabili per recipienti disponibili per piattaforme Solaris 2000

Per piattaforme Solaris 4000

	No. art.	Universale		A due livelli	
		18 x 30	36 x 24	18 x 30	
Supporto per recipiente singolo regolabile in altezza 11"x14"	75004104	-	_	_	
Supporto per recipiente a doppia regolazione in altezza 12"x14"	75004102	2	2	4	
Supporto per recipiente a doppia regolazione in altezza 18"x18"	75004103	1	2	2	
Supporto regolabile per recipiente singolo	75004101	8	12	16	

Tabella 25: Supporti regolabili per recipienti per piattaforme Solaris 4000

Per piattaforme Solaris 2000 I / 2000 R

	No. art.	Universale
		11 x 14
Supporto per recipiente singolo regolabile in altezza 11"x14"	75004104	1
Supporto per recipiente a doppia regolazione in altezza 12"x14"	75004102	-
Supporto per recipiente a doppia regolazione in altezza 18"x18"	75004103	-
Supporto regolabile per recipiente singolo	75004101	2

Tabella 26: Supporti regolabili per recipienti per piattaforme Solaris 2000 I / 2000 R

Per piattaforme Solaris 4000 I / 4000 R

	No. art.	Universale
		18 x 18
Supporto per recipiente singolo regolabile in altezza 11"x14"	75004104	-
Supporto per recipiente a doppia regolazione in altezza 12"x14"	75004102	1
Supporto per recipiente a doppia regolazione in altezza 18"x18"	75004103	1
Supporto regolabile per recipiente singolo	75004101	4

Tabella 27: Supporti regolabili per recipienti per piattaforme Solaris 4000 I / 4000 R

1. 2. 7. Supporti per recipienti quadrati

Per piattaforme Solaris 2000



	No. art.	Universale			A due livelli	
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Supporto per bottiglia quadrata da 125 ml	75004106	12	16	24	21	32
Supporto per bottiglia quadrata da 250 ml	75004107	9	16	22	18	32
Supporto per bottiglia quadrata da 500 ml	75004108	6	9	12	11	18
Supporto per bottiglia quadrata da 1000 ml	75004109	4	9	12	8	18

Tabella 28: Supporti per recipienti quadrati disponibili per piattaforme Solaris 2000

Per piattaforme Solaris 4000

	No. art.	Universale		A due livelli
		18 x 30	36 x 24	18 x 30
Supporto per bottiglia quadrata da 125 ml	75004106	28	54	56
Supporto per bottiglia quadrata da 250 ml	75004107	28	52	56
Supporto per bottiglia quadrata da 500 ml	75004108	15	24	30
Supporto per bottiglia quadrata da 1000 ml	75004109	15	24	28

Tabella 29: Supporti per recipienti quadrati disponibili per piattaforme Solaris 4000

Per piattaforme Solaris 2000 I / 2000 R

	No. art.	Universale
		11 x 14
Supporto per bottiglia quadrata da 125 ml	75004106	9
Supporto per bottiglia quadrata da 250 ml	75004107	7
Supporto per bottiglia quadrata da 500 ml	75004108	3
Supporto per bottiglia quadrata da 1000 ml	75004109	2

Tabella 30: Supporti per recipienti quadrati disponibili per piattaforme Solaris 2000 I / 2000 R

Per piattaforme Solaris 4000 I / 4000 R

	No. art.	Universale
		18 x 18
Supporto per bottiglia quadrata da 125 ml	75004106	16
Supporto per bottiglia quadrata da 250 ml	75004107	16
Supporto per bottiglia quadrata da 500 ml	75004108	9
Supporto per bottiglia quadrata da 1000 ml	75004109	9

Tabella 31: Supporti per recipienti quadrati disponibili per piattaforme Solaris 4000 I / 4000 R

1. 2. 8. Rack porta becher Nalgene

Per piattaforme Solaris 2000



	No. art.	Universale		A due livelli		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Dimensioni 11 x 14 / 12 x 14			Rack / posti			
Becher Nalgene da 30 ml, 36 posti	75004116	1/36	-	-	2/72	-
Becher Nalgene da 50 ml, 28 posti	75004129	1 / 28	-	-	2/56	-
Becher Nalgene da 100 ml, 16 posti	75004118	1 / 16	-	-	2/32	-
Becher Nalgene da 250 ml, 9 posti	75004119	1/9	-	-	2/18	-
Becher Nalgene da 400 ml, 8 posti	75004120	1/8	-	_	2/16	-
Becher Nalgene da 600 ml, 5 posti	75004121	1/5	-	-	2/10	-
Dimensioni 9 x 18			F	Rack / posti		
Becher Nalgene da 30 ml, 32 posti	75004110	-	2 / 64	2/64	-	4 / 128
Becher Nalgene da 50 ml, 24 posti	75004128	-	2 / 48	2/48	-	4 / 96
Becher Nalgene da 100 ml, 15 posti	75004112	-	2 / 30	2/30	-	4 / 60
Becher Nalgene da 250 ml, 8 posti	75004113	-	2/16	2/16	-	4 / 32
Becher Nalgene da 400 ml, 6 posti	75004114	-	2/12	2/12	-	4 / 24
Becher Nalgene da 600 ml, 4 posti	75004115	-	2/8	2/8	-	4 / 16

Tabella 32: Rack porta becher Nalgene disponibili per piattaforme Solaris 2000

Per piattaforme Solaris 4000

	No. art.	Universale		A due livelli	
		18 x 30	36 x 24	18 x 30	
Dimensioni 11 x 14 / 12 x 14			Rack / posti		
Becher Nalgene da 30 ml, 36 posti	75004116	-	-	-	
Becher Nalgene da 50 ml, 28 posti	75004129	-	-	-	
Becher Nalgene da 100 ml, 15 posti	75004118	-	-	-	
Becher Nalgene da 250 ml, 9 posti	75004119	-	-	-	
Becher Nalgene da 400 ml, 8 posti	75004120	-	-	-	
Becher Nalgene da 600 ml, 5 posti	75004121	-	-	-	
Dimensioni 9 x 18			Rack / posti		
Becher Nalgene da 30 ml, 32 posti	75004110	3 / 96	4 / 128	6 / 192	
Becher Nalgene da 50 ml, 24 posti	75004128	3/72	4 / 96	6/144	
Becher Nalgene da 100 ml, 15 posti	75004112	3 / 45	4 / 60	6 / 90	
Becher Nalgene da 250 ml, 8 posti	75004113	3/24	4 / 32	6 / 48	
Becher Nalgene da 400 ml, 6 posti	75004114	3 / 18	4 / 24	6/36	
Becher Nalgene da 600 ml, 4 posti	75004115	3/12	4 / 16	6 / 24	

Tabella 33: Rack porta becher Nalgene disponibili per piattaforme Solaris 4000
Per piattaforme Solaris 2000 I / 2000 R

	No. art.	Universale
		11 x 14
Dimensioni 11 x 14 / 12 x 14		Rack / posti
Becher Nalgene da 30 ml, 36 posti	75004116	1 / 36
Becher Nalgene da 50 ml, 28 posti	75004129	1 / 28
Becher Nalgene da 100 ml, 16 posti	75004118	1 / 16
Becher Nalgene da 250 ml, 9 posti	75004119	1/9
Becher Nalgene da 400 ml, 8 posti	75004120	1/8
Becher Nalgene da 600 ml, 5 posti	75004121	1/5
Dimensioni 9 x 18		Rack / posti
Becher Nalgene da 30 ml, 32 posti	75004110	-
Becher Nalgene da 50 ml, 24 posti	75004128	-
Becher Nalgene da 100 ml, 15 posti	75004112	-
Becher Nalgene da 250 ml, 8 posti	75004113	-
Becher Nalgene da 400 ml, 6 posti	75004114	-
Becher Nalgene da 600 ml, 4 posti	75004115	-

Tabella 34: Rack porta becher Nalgene disponibili per R piattaforme Solaris 2000 I / 2000 R

Per piattaforme Solaris 4000 I / 4000 R

	No. art.	Universale
		18 x 18
Dimensioni 11 x 14 / 12 x 14		Rack / posti
Becher Nalgene da 30 ml, 36 posti	75004116	-
Becher Nalgene da 50 ml, 28 posti	75004129	-
Becher Nalgene da 100 ml, 16 posti	75004118	-
Becher Nalgene da 250 ml, 9 posti	75004119	-
Becher Nalgene da 400 ml, 8 posti	75004120	-
Becher Nalgene da 600 ml, 5 posti	75004121	-
Dimensioni 9 x 18		Rack / posti
Becher Nalgene da 30 ml, 32 posti	75004110	2 / 64
Becher Nalgene da 50 ml, 24 posti	75004128	2 / 48
Becher Nalgene da 100 ml, 15 posti	75004112	2 / 30
Becher Nalgene da 250 ml, 8 posti	75004113	2/16
Becher Nalgene da 400 ml, 6 posti	75004114	2/12
Becher Nalgene da 600 ml, 4 posti	75004115	2/8

Tabella 35: Rack porta becher Nalgene disponibili per R piattaforme Solaris 4000 I / 4000 R

1. 2. 9. Supporto per imbuto separatore

Per piattaforme Solaris 2000



	No. art.	Universale			A due livelli		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18	
Supporto singolo per imbuto separatore, da 250 ml a 2000 ml	75004125	2	3	4	2	5	
Supporto per imbuto separatore, a montaggio verticale		4	4	6	-	-	

Tabella 36: Supporti per imbuto separatore disponibili per piattaforme Solaris 2000

Per piattaforme Solaris 4000

	No. art.	Unive	rsale	A due livelli	
		18 x 30	36 x 24	18 x 30	
Supporto singolo per imbuto separatore, da 250 ml a 2 000 ml	75004125	5	7	9	
Supporto per imbuto separatore, a montaggio verticale		8	12	-	

 Tabella 37:
 Supporti per imbuto separatore disponibili per piattaforme Solaris 4000

Per piattaforme Solaris 2000 I / 2000 R

	No. art.	Universale
Supporto singolo per imbuto separatore, da 250 ml a 2000 ml	75004125	-
Supporto per imbuto separatore, a montaggio verticale		-

Tabella 38: Supporti per imbuto separatore disponibili per piattaforme Solaris 2000 I / 2000 R

Per piattaforme Solaris 4000 I / 4000 R

	No. art.	Universale
		18 x 18
Supporto singolo per imbuto separatore, da 250 ml a 2000 ml	75004125	-
Supporto per imbuto separatore, a montaggio verticale		-

Tabella 39: Supporti per imbuto separatore disponibili per piattaforme Solaris 4000 I / 4000 R

1. 2. 10. Supporto multiuso

Per piattaforme Solaris 2000



	No. art.	Universale		A due livelli		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Supporto multiuso 11 x 14 / 12 x 14	75004123	1	-	-	2	-
Supporto multiuso 9 x 18	75004122	-	2	З	-	4

Tabella 40: Supporti multiuso disponibili per piattaforme Solaris 2000

Per piattaforme Solaris 4000

	No. art.	Unive	rsale	A due livelli	
		18 x 30	36 x 24	18 x 30	
Supporto multiuso 11 x 14 / 12 x 14	75004123	-	-	-	
Supporto multiuso 9 x 18	75004122	З	6	6	

Tabella 41: Supporti multiuso disponibili per piattaforme Solaris 4000

Per piattaforme Solaris 2000 I / 2000 R

	No. art.	Universale
		11 x 14
Supporto multiuso 11 x 14 / 12 x 14	75004123	1
Supporto multiuso 9 x 18	75004122	-

Tabella 42: Supporti multiuso disponibili per piattaforme Solaris 2000 I / 2000 R

Per piattaforme Solaris 4000 I / 4000 R

	No. art.	Universale
		18 x 18
Supporto multiuso 11 x 14 / 12 x 14	75004123	-
Supporto multiuso 9 x 18	75004122	2

Tabella 43: Supporti multiuso disponibili per piattaforme Solaris 4000 I / 4000 R

Accessori per supporto multiuso

	No. art.	Universale
Guamizioni O-ring di ricambio (supporto multiuso)	75004132	-

Tabella 44: Accessori disponibili per supporti multiuso

1. 2. 11. Tappetini antiscivolo

Informazioni generali sulla scelta dei tappetini antiscivolo:

- I tappetini con una grandezza di 170 x 280 mm sono ottimi per piattaforme da 11 x 14 e 12 x 14.
- I tappetini con una grandezza di 200 x 200 mm sono ottimi per piattaforme più grandi di dimensione 18 x 18, 18 x 24, 30 x 18 e 36 x 24.
- Tutti i tappetini possono essere tagliati a qualsiasi misura desiderata.

Per piattaforme Solaris 2000

	No. art.	Universale		A due livelli, universale		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
Tappetino 200 x 200 mm ad alta adesione, confezione da 2	75004126	1	4	6	2	8
Tappetino 280 x 170 mm ad alta adesione, confezione da 2	75004127	2	2	4	4	4
Tappetino 200 x 200 mm a bassa adesione, confezione da 2	75004111	1	4	6	2	8
Tappetino 280 x 170 mm a bassa adesione, confezione da 2	75004117	2	2	4	4	4

I tappetini sono venduti in confezioni da 2. Il numero di tappetini che trovano posto su una piattaforma non fa riferimento alla grandezza della confezione ma al singolo tappetino.

Tabella 45: Tappetini antiscivolo disponibili per piattaforme Solaris 2000

Per piattaforme Solaris 4000

	No. art.	Unive	rsale	A due livelli
		18 x 30	36 x 24	18 x 30
Tappetino 200 x 200 mm ad alta adesione, confezione da 2	75004126	6	12	12
Tappetino 280 x 170 mm ad alta adesione, confezione da 2	75004127	6	9	12
Tappetino 200 x 200 mm a bassa adesione, confezione da 2	75004111	6	12	12
Tappetino 280 x 170 mm a bassa adesione, confezione da 2	75004117	6	9	12

I tappetini sono venduti in confezioni da 2. Il numero di tappetini che trovano posto su una piattaforma non fa riferimento alla grandezza della confezione ma al singolo tappetino.

Tabella 46: Tappetini antiscivolo disponibili per piattaforme Solaris 4000

Per piattaforme Solaris 2000 I / 2000 R

	No. art.	Universale
Tappetino 200 x 200 mm ad alta adesione, confezione da 2	75004126	1
Tappetino 280 x 170 mm ad alta adesione, confezione da 2	75004127	2
Tappetino 200 x 200 mm a bassa adesione, confezione da 2	75004111	1
Tappetino 280 x 170 mm a bassa adesione, confezione da 2	75004117	2

I tappetini sono venduti in confezioni da 2. Il numero di tappetini che trovano posto su una piattaforma non fa riferimento alla grandezza della confezione ma al singolo tappetino.

Tabella 47: Tappetini antiscivolo disponibili per piattaforme Solaris 2000 I / 2000 R

Per piattaforme Solaris 4000 I / 4000 R

	No. art.	Universale
		18 x 18
Tappetino 200 x 200 mm ad alta adesione, confezione da 2	75004126	4
Tappetino 280 x 170 mm ad alta adesione, confezione da 2	75004127	2
Tappetino 200 x 200 mm a bassa adesione, confezione da 2	75004111	4
Tappetino 280 x 170 mm a bassa adesione, confezione da 2	75004117	2

I tappetini sono venduti in confezioni da 2. Il numero di tappetini che trovano posto su una piattaforma non fa riferimento alla grandezza della confezione ma al singolo tappetino.

Tabella 48: Tappetini antiscivolo disponibili per piattaforme Solaris 4000 I / 4000 R

1. 2. 12. Accessori generali

I. 2. 12. Accessori generali		
	No. art.	
Rullo applicatore per tappetini antiscivolo	75004124	
Cacciavite per accessori	75004131	
Collettore del gas (Solaris 2000 I / 2000 R)	SK2000-8GM	
Collettore del gas (Solaris 4000 I / 4000 R)	SK4000-8GM	
Kit di ricambi per supporto (viti)	SK0010	
Kit di ricambi piattaforma per Solaris 2000, 2000 VR e 4000 VR (viti per piattaforma, attrezzo)	SK0100	
Kit di ricambi piattaforma per SK4000 (viti per piattaforma, attrezzo)	SK0101	

Tabella 49: Accessori generali disponibili

1.3. Direttive e norme

Regione	Direttiva	Norme
Europa	2006/42/CE	EN 61010-1
	Direttiva Macchine	IEC 61010-2-051
	2011/65/EU RoHS	EN 61326-1 Classe B
	Restrizione dell'uso di	EN ISO 14971
	determinate sostanze pericolose	EN ISO 9001
	ed elettroniche	Solo per Solaris 2000 I / 2000 R /
	<u>Scopi protettivi:</u>	4000 I / 4000 R: IEC 61010-2-010
	2014/35/EU Bassa tensione	
	2014/30/CE Compatibilità	
	elettromagnetica (EMC)	
Nord America		ANSI/UL 61010-1
		IEC 61010-2-051
		IEC 61326-1 Classe B
		CFR 47 FCC 15 EMC
		EN ISO 14971
		EN ISO 9001
		Solo per Solaris 2000 I / 2000 R /
		4000 I / 4000 R: IEC 61010-2-010
Giappone		IEC 61010-1
		IEC 61010-2-051
		IEC 61326-1 Classe B
		EN ISO 14971
		EN ISO 9001
		Solo per Solaris 2000 I / 2000 R / 4000 I / 4000 R: IEC 61010-2-010

Tabella 50:Direttive e norme

AVVISO: Questo dispositivo è stato sottoposto a test ed è risultato conforme ai limiti previsti per i dispositivi digitali di Classe A ai sensi della Parte 15 della normativa FCC. Questi limiti sono stati stabiliti al fine di fornire una protezione adeguata da interferenze dannose quando il dispositivo viene utilizzato in un ambiente commerciale. Questo dispositivo genera, usa e può irradiare energia a radiofrequenza e, se non istallato e usato in conformità con le procedure descritte nel manuale d'uso, può causare interferenze dannose alle comunicazioni radio. È probabile che l'utilizzo di questo dispositivo in un'area residenziale provochi interferenze dannose. In tal caso, l'utente sarà tenuto a correggere le interferenze a proprie spese.

1.4. Informazioni generali sul prodotto

1.4.1. Solaris 2000



① Piastra di montaggio piattaforma; ② Interfaccia utente grafica (GUI); ③ Interruttore generale; ④ Collegamento alla rete; ⑤ Fusibile; ⑥ Porta Ethernet; ⑦ Porta USB

Figura 1: Solaris 2000 - Vista d'insieme

1.4.2. Solaris 4000



① Piastra di montaggio piattaforma; ② Interfaccia utente grafica (GUI); ③ Interruttore generale; ④ Collegamento alla rete; ⑤ Fusibile (ripristinabile); ⑥ Porta Ethernet; ⑦ Porta USB

Figura 2: Solaris 4000 - Vista d'insieme

1.4.3. Solaris 2000 I / 2000 R



① Piastra di montaggio piattaforma; ② Interfaccia utente grafica (GUI); ③ Interruttore generale; ④ Foro passacavo; ⑤ Collegamento alla rete; ⑥ Fusibile (ripristinabile); ⑦ Porta Ethernet; ⑧ Porta USB; ⑨ Cappa

Figura 3: Vista d'insieme Solaris 2000 I / 2000 R

1.4.4. Solaris 4000 I / 4000 R



Piastra di montaggio piattaforma;
 Interfaccia utente grafica (GUI);
 Interruttore generale;
 Foro passacavo;
 Collegamento alla rete;
 Fusibile (ripristinabile);
 Porta Ethernet;
 Porta USB;
 Cappa

Figura 4: Vista d'insieme Solaris 4000 I / 4000 R

1.4.5. Raccordi

Agitatori aperti



① Collegamento alla rete; ② Fusibile; ③ Porta Ethernet

Figura 5: Vista posteriore (a sinistra: Solaris 4000, a destra: Solaris 2000)

Agitatori-incubatori (riscaldati) e agitatori refrigerati (raffreddati e riscaldati)



① Collegamento alla rete; ② Fusibile; ③ Porta Ethernet; ④ Tubo di drenaggio
 Figura 6: Vista posteriore (a sinistra: Solaris 2000 I / 2000 R, a destra: Solaris 4000 I / 4000 R)

Collegamento alla rete

L'agitatore richiede una fonte di alimentazione elettrica conforme alle proprie specifiche. I cavi di alimentazione sono in dotazione.

No. art.	Agitatore	Specifiche
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

Figura 7: Specifiche per l'alimentazione elettrica degli agitatori

AVVERTENZA Danno causato da un'alimentazione elettrica o da una spina elettrica non appropriate. Assicurarsi che l'agitatore sia collegato a una presa correttamente messa a terra secondo le norme di sicurezza elettrica vigenti. Non usare l'agitatore con un cavo di alimentazione difettoso o inadeguato.

La spina di rete deve essere sempre liberamente accessibile.

Per evitare il rischio di scosse elettriche, assicurarsi che la superficie intorno all'agitatore sia asciutta. In caso di versamento accidentale o di spruzzi di liquidi, sconnettere l'agitatore dall'alimentazione elettrica, rimuovere i liquidi traboccati e eliminare qualsiasi rischio biologico o di salute prima di continuare.

Sconnettere l'agitatore dalla rete di alimentazione quando non viene utilizzato.

Fusibile

Il fusibile scattato a seguito di condizioni di sovracorrente può essere ripristinato. L'agitatore non sarà operativo finché il fusibile non sia stato riportato nella sua posizione corretta.

Ethernet

L'agitatore dispone di una porta Ethernet RJ45 che può essere utilizzata per il collegamento alla Local Area Network (LAN). Usare solo apparecchiature conformi alla norma IEC 60950-1 con la porta Ethernet RJ45. La porta Ethernet RJ45 permette di caricare futuri aggiornamenti del software.

Fori passacavo

Gli agitatori a termoregolazione dispongono di due aperture laterali denominate fori passacavo. Questi fori passacavo sono solitamente chiusi con dei tappi. Il tappo può essere rimosso per introdurre strumenti aggiuntivi, come ad esempio un ulteriore sensore di temperatura, oppure per passare cavi o tubi, come i tubi del collettore del gas.

<u>USB</u>

L'agitatore è dotato di due porte USB-A 2.0 che possono essere utilizzate con una chiavetta USB comunemente disponibile in commercio. Usare solo apparecchiature conformi alla norma IEC 60950-1 con le porte USB.

Protezione interna da sovratemperatura

Gli agitatori a termoregolazione dispongono di un termostato con ripristino manuale installato accanto agli elementi riscaldanti. Questo dispositivo termico spegne i riscaldatori per evitare temperature eccessivamente elevate all'interno della camera dell'agitatore in caso di un guasto ai ventilatori di circolazione aria.

Il termostato a reset manuale deve essere resettato manualmente per ripristinare il funzionamento dell'agitatore. Ciò può essere effettuato solo dal servizio di assistenza di Thermo Fisher Scientific.

2. Trasporto e installazione



L'operatore stesso è responsabile della corretta installazione dell'agitatore.

AVVISO

Si consiglia di controllare il cartone d'imballaggio al momento della consegna. Prima di spacchettare il dispositivo controllare l'imballaggio per individuare eventuali danni causati durante il trasporto. Se sono riscontrati danni, lo spedizioniere deve specificare il danno e firmare la copia della bolletta di consegna per confermare il danno.

Aprire con cautela il cartone e assicurarsi che siano presenti tutti i componenti («Tabella 51: Dotazione di fornitura») prima di smaltire il materiale d'imballaggio. Se dopo il disimballo sono riscontrati danni, informare lo spedizioniere e richiedere un'ispezione del danno.

Importante: Se entro pochi giorni dalla consegna del prodotto non avviene la richiesta di ispezione del danno, lo spedizioniere sarà esonerato da qualsiasi responsabilità per il danno. Deve essere chiesta un'ispezione del danno.

2.1. Apertura dell'imballo

Usare la distinta della merce spedita per verificare che l'unità sia stata consegnata completa. Non smaltire il materiale d'imballaggio prima di aver individuato tutti i componenti.

Dotazione di fornitura

Pos.	Quantità
Agitatore	1
Cavo di alimentazione	1
Piattaforma universale	1*
Viti per piattaforma	
- Solaris 4000	4*
- altri modelli	3*
Manuale inglese in forma cartacea	1
Manuali su chiavetta USB	1
Attrezzo di fissaggio (chiave a T)	1
* Gli agitatori 2000 I, 2000 R, 4000 I e 4000 R preinstallata nell'agitatore. Né la piattaforma né come componenti separati.	hanno la piattaforma già le viti sono disponibili

Tabella 51: Dotazione di fornitura

Se mancano dei componenti, rivolgersi a Thermo Fisher Scientific.

2.2. Luogo di installazione

ATTENZIONE La protezione può essere pregiudicata a seguito di una stabilità ridotta delle materie plastiche esposte ai raggi ultravioletti. Non esporre l'agitatore e gli accessori di plastica ai raggi diretti del sole o ad altri raggi ultravioletti.

Posizionare l'agitatore su un tavolo o banco piano in grado di reggere il peso dell'agitatore, compresi gli accessori e campioni, durante il funzionamento. Collocare l'agitatore in prossimità di una presa elettrica che soddisfi i requisiti riportati sulla targhetta. Lasciare uno spazio sufficiente intorno al dispositivo per garantire la libera circolazione dell'aria, uno spazio sufficiente per gli accessori e un maggiore comfort per l'utente.

Attenersi ai seguenti requisiti per l'installazione del dispositivo:

- Per assicurare una distanza sufficiente verso parti mobili ed evitare lesioni da schiacciamento o danni ad attrezzature adiacenti, l'agitatore aperto richiede uno spazio libero di 8 cm (3") su tutti i lati della piattaforma. La piattaforma dell'agitatore aperto può sporgere oltre il corpo.
- Gli agitatori causano vibrazioni. Non conservare dispositivi sensibili oppure sostanze o oggetti pericolosi nella zona di sicurezza.



Figura 8: Collocazione dell'agitatore aperto con uno spazio libero di 8 cm (3")

Per garantire una ventilazione accurata e ottime prestazioni degli agitatori a termoregolazione, è necessario rispettare le seguenti distanze:

- Gli agitatori-incubatori (riscaldati) Solaris 2000 I e 4000 I richiedono uno spazio libero di 8 cm (3") su tutti i quattro lati del corpo, come illustrato sul lato sinistro della Figura 9.
- Gli agitatori refrigerati (raffreddati e riscaldati) Solaris 2000 R e 4000 R possono essere usati con uno spazio libero di 8 cm (3") su tutti i quattro lati del corpo se sopra la cappa vi è uno **spazio libero** come illustrato sul lato sinistro della Figura 9.

 Se lo spazio libero sopra l'agitatore refrigerato (raffreddato e riscaldato) Solaris 2000 R e 4000 R è ostacolato, lo spazio libero intorno al corpo deve essere aumentato a 30 cm (12") su tutti i quattro lati, come illustrato sul lato destro della Figura 9.



Figura 9: Spazi liberi per la collocazione degli agitatori-incubatori (riscaldati) (a sinistra) e per gli agitatori refrigerati (raffreddati e riscaldati) (a destra); per gli agitatori refrigerati (raffreddati e riscaldati) con uno spazio libero sufficiente verso l'alto vale la figura a sinistra.

AVVERTENZA Durante l'agitazione eventuali sostanze pericolose devono essere tenute fuori da questa zona.

Il piano di appoggio deve:

- » essere pulito,
- » essere stabile, solido, rigido ed esente da vibrazioni,
- » permettere il posizionamento orizzontale dell'agitatore,
- » sopportare il peso dell'agitatore.
- L'agitatore deve essere utilizzato solo in ambienti interni.
- L'agitatore non deve essere esposto a fonti di calore né tanto meno a forte luce solare.
- Il luogo di installazione deve essere sempre ben aerato.
- La spina di rete deve essere sempre liberamente accessibile.

2.3. Trasporto

ATTENZIONE Lesioni causate dalla caduta dell'agitatore dall'alto. Sollevare sempre l'agitatore da entrambi i lati. Non sollevare mai l'agitatore, afferrandolo per il pannello frontale o per la piattaforma installata. Non sollevare mai un agitatore a termoregolazione afferandolo per la maniglia della cappa trasparente.

ATTENZIONE L'uso sicuro dell'agitatore viene pregiudicato dalla condensazione di acqua all'interno del corpo. Ciò accade quando l'agitatore viene portato da un posto più fresco in uno più caldo con aria più umida, ad esempio dall'area di immagazzinamento o dal camion di trasporto nella sala laboratorio. Lasciare asciugare l'apparecchiatura per almeno 2 ore prima dell'utilizzo.

AVVISO Con gli agitatori aperti rimuovere sempre la piattaforma, il carico e tutti gli accessori prima di spostare l'agitatore. Se la piattaforma non viene rimossa, sussiste il pericolo di danneggiare la piastra di montaggio o il meccanismo di agitazione. Con gli agitatori a termoregolazione è sufficiente rimuovere il carico dalla piattaforma.

AVVISO Il dispositivo può essere danneggiato da urti. Trasportare l'agitatore in posizione verticale e se possibile con l'imballaggio originale.

Manipolazione dell'agitatore

Durante la manipolazione dell'agitatore fare attenzione a quanto segue:

- Sollevare l'agitatore dai due lati, ma mai da davanti e dietro
- Impiegare il numero di persone necessarie in considerazione del peso dell'agitatore (per maggiori informazioni consultare «Specifiche tecniche» a pagina 12), comunque mai meno di due persone





Figura 10: Come trasportare un agitatore

AVVERTENZA Sollevare sempre l'agitatore da entrambi i lati. Non sollevare mai l'agitatore afferrandolo sul lato frontale e posteriore. Gli agitatori sono pesanti (per maggiori informazioni consultare «Specifiche tecniche» a pagina 12). Sollevare e trasportare l'agitatore con almeno due persone.

ATTENZIONE Per trasportare un agitatore a termoregolazione non afferrarlo mai per la maniglia della cappa trasparente. La maniglia potrebbe rompersi e l'agitatore potrebbe cadere a terra, provocando gravi lesioni alle persone.

2.4. Livellamento

L'agitatore stesso non dispone di una possibilità di livellamento. Per un'installazione accurata dell'agitatore serve un piano di appoggio livellato.

AVVISO Non posizionare spessori o altri oggetti piatti tra i piedi dell'agitatore e il piano di appoggio per livellare l'agitatore.

2.5. Collegamento alla rete

L'agitatore richiede una fonte di alimentazione elettrica conforme alle proprie specifiche. I cavi di alimentazione sono in dotazione. Per maggiori informazioni consultare la Figura 7 a pagina 48.

AVVERTENZA Danno causato da un'alimentazione elettrica o da una spina non corrette. Assicurarsi che l'agitatore sia collegato sempre solo a prese correttamente messe a terra. Non usare l'agitatore con un cavo di alimentazione difettoso o inadeguato.

AVVISO Radiazioni elettromagnetiche possono causare interferenze sul display. Questa circostanza non danneggia il dispositivo e non limita o altera la sua funzionalità. Per evitare interferenze provocate da radiazioni elettromagnetiche si consiglia di non portare dispositivi mobili, come i cellulari, nelle immediate vicinanze del dispositivo. Non fare funzionare il dispositivo insieme ad altri dispositivi ad alta potenza elettrica sullo stesso circuito. Non usare più dispositivi collegati alla stessa presa multipla.

Per il collegamento dell'agitatore alla rete di alimentazione seguire la seguente procedura:

- 1. Disinserire l'interruttore generale che si trova sul lato destro dell'apparecchiatura.
- 2. Assicurarsi che le specifiche del cavo corrispondano allo standard vigente nel proprio Paese.
- 3. Assicurarsi che la tensione e la frequenza corrispondano ai dati riportati sulla targhetta.

La spina di rete deve essere sempre liberamente accessibile.

Per evitare il rischio di scosse elettriche, assicurarsi che la superficie intorno all'agitatore sia asciutta. In caso di versamento accidentale o di spruzzi di liquidi, sconnettere l'agitatore dall'alimentazione elettrica ed eliminare i liquidi traboccati prima di continuare.

Sconnettere l'agitatore dalla rete di alimentazione quando non viene utilizzato.

2.6. Configurazione iniziale

Eseguire le seguenti operazioni prima di utilizzare il dispositivo:

1. Accendere il dispositivo. Viene visualizzato il logo di Thermo Scientific. Toccare il pulsante **Start Setup**.



Figura 11: Richiesta di configurazione iniziale

2. Selezionare la lingua desiderata sulla schermata Lingua. Toccare il pulsante Avanti.



Figura 12: Configurazione iniziale - Lingua

 Se desiderato, è possibile inserire il nome del dispositivo nella finestra di dialogo Nome dell'unità. Toccare il pulsante Avanti.

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

Figura 13: Configurazione iniziale - Nome dell'unità

 Se desiderato, è possibile inserire la città e il paese nel campo di testo. Dopo aver digitato tre caratteri, nel campo di testo per la città e il paese appare un elenco con suggerimenti. Se desiderato, scegliere una voce e toccare il pulsante Avanti.

Set Region			
	Enter the city and c:	untry closest to your time zone	
City, Country:	Tap to ente	r	
B	ack	Next	

Figura 14: Configurazione iniziale - Regione

5. Selezionare il formato data. Toccare il pulsante Avanti.

Set Date Disp	lay Format		
O YYYY/MMM/DD			
Back	Next		

Figura 15: Configurazione iniziale - Imposta formato di visualizzazione data

6. Ruotare i singoli quadranti della rotella di selezione per selezionare la data attuale. Toccare il pulsante **Avanti.**



Figura 16: Configurazione iniziale - Imposta Data

7. Selezionare il formato orario e ruotare i singoli quadranti della rotella di selezione per impostare l'ora attuale. Toccare il pulsante **Avanti**.



Figura 17: Configurazione iniziale - Imposta Ora

 Solo per agitatori-incubatori (riscaldati) e agitatori refrigerati (raffreddati e riscaldati): Selezionare l'unità di misura desiderata per la temperatura. Toccare il pulsante Avanti.

Temperat	ture Unite		
rempera	ture onito		
	*0		
•	-0		
O *E			
0			
Back	Next		

Figura 18: Configurazione iniziale - Selezione dell'unità di misura per la temperatura

 Solo per agitatori-incubatori (riscaldati) e agitatori refrigerati (raffreddati e riscaldati): Cliccare sulla freccia blu per impostare l'allarme per temperatura alta e i valori limite per un arresto d'emergenza. Toccare il pulsante Avanti.

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

Figura 19: Configurazione iniziale - Impostazione dei valori limite per l'allarme di temperatura alta e bassa

10. Ripetere la procedura per i valori limite per l'allarme di temperatura bassa. Toccare **Avanti** per continuare.

- 11. Se si desidera continuare senza impostare un codice di accesso da amministratore, lasciare **Open Mode** (senza controllo dell'accesso) e toccare
 - Avanti per continuare.

Se si desidera impostare un codice di accesso da amministratore, toccare **Secure Mode** (con controllo dell'accesso).

•	Open Mode			
	Select to allow full access to all functionality for all users.			
0	Secure Mode			
	Select to require users to enter the admin passcode to make			
	(All users still able to use programs and set new run parameter	ers.		
		-		

Figura 20: Configurazione iniziale - Selezione della modalità di accesso

12. Nella finestra di richiesta del codice di accesso che appare, inserire un codice di accesso da amministratore con la tastiera e toccare **Avanti**.



Figura 21: Configurazione iniziale - Selezione della modalità di accesso

- 13. Nella successiva finestra di richiesta del codice di accesso, inserire il codice da amministratore.
- 14. Nella terza richiesta di inserimento del codice di accesso inserire ancora una volta il codice da amministratore per la conferma.



Figura 22: Configurazione iniziale - Inserimento e conferma del nuovo codice di accesso da amministratore

15. Sulla schermata di conferma che appare, toccare Avanti per continuare.

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				

Figura 23: Configurazione iniziale - Conferma della modifica al codice di accesso da amministratore

16. Viene visualizzata una schermata con istruzioni generali per l'installazione fisica dell'agitatore. Toccare il pulsante **Avanti**.

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				

Figura 24: Configurazione iniziale - Istruzioni per l'installazione

17. Viene visualizzata la schermata di configurazione completata. Toccare **Avanti** per concludere la procedura.



Figura 25: Configurazione iniziale - Calibrazione o fine della configurazione

AVVISO Sugli agitatori-incubatori (riscaldati) e sugli agitatori refrigerati (raffreddati e riscaldati) si può toccare **Calibra** e avviare la procedura descritta al capitolo «4. 6. Calibrazione di temperatura» a pagina 127.

2.7. Stoccaggio



Se l'agitatore e gli accessori sono messi fuori servizio, l'intero sistema deve essere pulito e, se necessario, disinfettato o decontaminato. Non lasciare l'agitatore e gli accessori in uno

ATTENZIONE stato di contaminazione indefinito. In caso di dubbi riguardanti la procedura da seguire, contattare il servizio di assistenza Thermo Fisher Scientific («Pulizia» a pagina 124, «Disinfezione» a pagina 125 e «Decontaminazione» a pagina 125).

- Prima di immagazzinare l'agitatore e gli accessori, questi devono essere puliti e, all'occorrenza, disinfettati e decontaminati.
- L'agitatore e gli accessori devono essere completamente asciutti prima dell'immagazzinamento.
- Conservare l'agitatore in un luogo pulito e privo di polvere.
- Appoggiare l'agitatore sui propri piedi.
- Non conservare l'agitatore in un luogo con esposizione diretta ai raggi del sole.

2.8. Spedizione



ATTENZIONE

Prima di spedire l'agitatore e gli accessori, l'intero sistema deve essere pulito e, se necessario, disinfettato o decontaminato. Non lasciare l'agitatore e gli accessori in uno stato di contaminazione indefinito. In caso di dubbi riguardanti la procedura da seguire, contattare il servizio di assistenza Thermo Fisher Scientific («Pulizia» a pagina 124, «Disinfezione» a pagina 125 e «Decontaminazione» a pagina 125).

Prima di spedire l'agitatore:

- L'agitatore deve essere pulito e decontaminato.
- La decontaminazione deve essere confermata con un certificato di decontaminazione. Un certificato di decontaminazione può essere richiesto al servizio di assistenza ai clienti di Thermo Fisher Scientific.

3. Funzionamento

3.1. Accensione/spegnimento

Premere l'interruttore generale sul lato destro del dispositivo per accendere (I) o spegnere (0) l'agitatore.

Durante la procedura di avvio sul touch-screen appare il logo di Thermo Scientific.

Quindi, il touch-screen visualizza l'attuale stato dell'agitatore.

3.2. Interfaccia utente grafica

La schermata principale (Home) è la schermata standard per l'interfaccia utente grafica (GUI) dell'agitatore. Rappresenta il punto di partenza per l'uso dell'agitatore. La schermata principale permette:

- di impostare i parametri operativi base, come ad esempio la velocità, la durata e la temperatura di agitazione (solo per agitatori a termoregolazione)
- di avviare e arrestare l'agitatore
- di visualizzare le informazioni di stato e di gestire gli allarmi e gli avvertimenti
- di passare ad altre schermate con informazioni di stato e opzioni di configurazione

Il contenuto della schermata differisce leggermente tra i modelli di agitatore aperto e quelli a termoregolazione.

La schermata principale per l'agitatore aperto si presenta come l'esempio illustrato nella Figura 26.



Figura 26: Schermata principale dell'interfaccia utente per agitatori aperti

La schermata principale per l'agitatore a termoregolazione dispone inoltre di opzioni di regolazione della temperatura, come illustrato nella Figura 27.



Figura 27: Schermata principale dell'interfaccia utente per agitatori a termoregolazione

3. 2. 1. Impostazione dei parametri operativi base per

l'agitatore

Sulla schermata principale basta un solo tocco per accedere alle schermate di configurazione dei parametri operativi più importanti dell'agitatore.

Impostazione della velocità

1. Toccare il campo Velocità attuale (© nella Figura 26 o ⑦ nella Figura 27) per accedere alla schermata del Valore nominale di velocità illustrato nella Figura 28.



Figura 28: Valore nominale di velocità

- Toccare le frecce sopra o sotto ogni quadrante della rotella di selezione per impostare la velocità desiderata. In alternativa, è anche possibile modificare la velocità nominale desiderata, ruotando il singolo quadrante della rotella di selezione.
- 3. Toccare il pulsante **Salva**.

AVVISO Se appare una finestra popup con l'informazione che la velocità nominale è fuori campo, il valore impostato non rientra nel campo di velocità supportato dall'agitatore, come specificato al capitolo «Dati tecnici» a pagina 12. Correggere l'impostazione e continuare.

4. Confermare la finestra di dialogo **Le impostazioni sono state salvate** che appare per ritornare alla schermata principale.

Calcolatore orbita: Calcolo approssimativo dalle impostazioni precedenti del dispositivo

Nella finestra di dialogo del valore nominale di velocità si può toccare anche il pulsante **Calcola velocità** per accedere al **Calcolatore orbita**: Il calcolatore orbita è utile quando si passa da un agitatore di struttura meccanica diversa a uno degli agitatori della serie Solaris descritti nel presente manuale. Questo calcolatore usa la differenza nella dimensione dell'orbita per dare una stima approssimativa della velocità nominale richiesta per ottenere risultati simili.

AVVISO Non usare direttamente i risultati ottenuti con il calcolatore orbita per la produzione dei campioni, ma effettuare una serie di prove per confermare i valori. In base agli sforzi di taglio presenti, le cellule possono crescere in maniera più rapida o più lenta, esprimere proteine diverse o essere danneggiate (o morire) dato che le condizioni fisiche cambiano con lo spostamento in un'orbita diversa.

1. Selezionare l'unità di misura in millimetri o pollici.



Figura 29: Calcolatore orbita

 Selezionare la dimensione dell'orbita precedente. Toccare il pulsante Calcola. Durante il calcolo della velocità viene visualizzata la seguente finestra.



Figura 30: Orbita calcolata

- 3. Toccare il pulsante **OK**.
- 4. Può essere calcolata la velocità consigliata per una dimensione personalizzata dell'orbita. Per effettuare il calcolo per un'orbita personalizzata, selezionare **Altro**.
- 5. Inserire la dimensione dell'orbita.



Figura 31: Calcolo dell'orbita di dimensione personalizzata

6. Toccare il pulsante Calcola.

Appare la finestra Velocità calcolata, visualizzando la velocità per la dimensione inserita.

Impostazione della durata del ciclo

L'agitatore può essere usato in modalità continua o temporizzata. In modalità continua l'agitatore può essere fermato manualmente a propria discrezione. In modalità temporizzata l'azionamento dell'agitatore si ferma automaticamente dopo che è trascorso il tempo impostato. Il temporizzatore può essere visualizzato come segue:

- Tempo trascorso: tempo trascorso, cioè per quanto tempo l'agitatore sta già funzionando dal momento in cui è stato premuto il pulsante di avvio, oppure
- Tempo rimanente: tempo residuo, cioè per quanto tempo l'agitatore continuerà a funzionare fino al trascorrere del tempo impostato.

- 1. Toccare il campo del **Tempo** (5) nella Figura 26 e Figura 27) per accedere alla schermata Modalità Tempo illustrata nella Figura 32.
- 2. Scegliere **Temporizzato**.



Figura 32: Selezione della Modalità Tempo

- 3. Inserire le ore e i minuti nelle caselle Imposta Tempo.
- 4. Toccare il pulsante Salva.
- 5. Per ritornare alla schermata principale confermare la finestra di dialogo Le impostazioni sono state salvate che appare.

Impostazione della temperatura

Sui modelli a termoregolazione, il campo **Temp nominale** permette di impostare, in qualsiasi momento, una temperatura per la rispettiva applicazione.

- Toccare il campo **Temp nominale** (parte inferiore della posizione

 nella Figura
 per accedere alla schermata **Temp nominale** illustrata nella Figura 33.
- Toccare le frecce sopra o sotto ogni quadrante della rotella di selezione per preimpostare la temperatura standard che l'agitatore deve mantenere. In alternativa, è anche possibile modificare la temperatura desiderata, ruotando il singolo quadrante della rotella di selezione.



Figura 33: Impostazione del valore nominale per la temperatura: schermata con l'indicazione della calibrazione

3. Toccare il pulsante **Salva**.

AVVISO Se appare una finestra popup con l'informazione che la temperatura nominale è fuori campo, il valore impostato non rientra nel campo di temperatura supportato dall'agitatore, come specificato al capitolo «Dati tecnici» a pagina 12. Correggere l'impostazione e continuare.

AVVISO Se appare una finestra popup con l'informazione che la temperatura selezionata potrebbe essere fuori campo a seguito dell'attuale temperatura ambiente, la temperatura ambiente non rientra nel campo di temperatura operativa dell'agitatore, come specificato al capitolo «Dati tecnici» a pagina 12. Toccare **OK** per confermare di aver letto il messaggio e continuare oppure selezionare una temperatura diversa.

AVVISO La schermata **Temp nominale** visualizza la temperatura calibrata e l'offset impostato durante la calibrazione, come illustrato nella parte in basso a sinistra della Figura 33. La procedura di calibrazione è descritta in «Calibrazione di temperatura» a pagina 127.

 Per ritornare alla schermata principale confermare la finestra di dialogo Le impostazioni sono state salvate che appare.

AVVISO Gli agitatori a termoregolazione hanno un comportamento specifico al raggiungere della temperatura impostata, chiamato oscillazione di temperatura. L'oscillazione di temperatura vuol dire che inizialmente la temperatura supera leggermente (o rimane leggermente sotto) la temperatura impostata per la camera e successivamente si avvicina alla temperatura impostata. Durante l'avvicinamento alla temperatura impostata il display non visualizza questo comportamento di oscillazione. Visualizza, invece, l'aumento (o la riduzione) della temperatura nella camera fino a raggiungere la temperatura impostata.

Avvio e arresto dell'agitatore

1. Per avviare l'agitatore toccare il pulsante Avvio.

Il pulsante Start cambia in Stop.



Figura 34: Pulsante di avvio e di arresto

2. Per arrestare l'agitatore toccare il pulsante Arresto.

3. 2. 2. Stato

Se l'agitatore si trova stato operativo regolare, sul touch-screen viene visualizzato il



simbolo di un cuore verde nell'area dello stato operativo e delle statistiche (7) nella Figura 26 e (8) nella Figura 27). Toccando il simbolo del cuore verde, si apre la schermata dello stato operativo. La schermata dello stato fornisce

informazioni statistiche sul funzionamento dell'agitatore, comprese le ore di agitazione e le ore operative totali nonché dell'ultima sessione. Gli agitatori a termoregolazione visualizzano, inoltre, le ore dei processi di raffreddamento e riscaldamento.

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	60brs	Hours Shaking	232hrs	Hours Shaking	60hrs
Haura Dawarad	2026		Clhus	Hours Powered	332hrs	Hours Powered	61hrs
Hours Powered	332nrs	Hours Powered	ornrs	Hrs tempering	168hrs	Hrs tempering	61hrs
			()				(i)

Figura 35: Dati statistici dell'agitatore: agitatore aperto (a sinistra) e agitatore a termoregolazione (a destra)

Avvertimento

Se viene emesso un avvertimento, sul display touch-screen appare una barra gialla con testo scorrevole in alto sullo schermo attuale. Inoltre, viene emesso un segnale acustico d'allarme. La barra gialla di avvertimento con testo

scorrevole scompare dopo aver fatto scorrere due volte il testo. Solo il triangolo giallo indica la presenza di uno o più avvertimenti per l'agitatore. Il simbolo della campana ha un cerchio celeste con bordo bianco che riporta il numero di avvertimenti attivi. Toccando il simbolo del triangolo nell'area di informazione e di stato operativo (⑦ nella Figura 26 e ⑧ nella Figura 27) si apre una schermata con un elenco di tutti gli avvertimenti attualmente attivi. L'ultimo avvertimento viene visualizzato con il testo completo per permettere all'operatore di vedere tutti i dettagli, come illustrato nella Figura 36. Si ha la possibilità di scorrere l'elenco e di toccare qualsiasi voce dell'elenco per estendere il testo e visualizzare maggiori dettagli.



Figura 36: Lista degli avvertimenti

Si ha la possibilità di selezionare l'avvertimento attivo, spuntando la casella accanto alla voce di avvertimento. Toccando il pulsante **Conferma** il sistema cerca di eliminare il messaggio di avvertimento dalla lista. Dopo l'eliminazione di tutti gli avvertimenti l'icona dello stato si ritrasforma nell'originario simbolo del cuore verde.

<u>Allarme</u>

Se viene emesso un allarme, l'agitatore si ferma subito per evitare danni ai campioni



e/o al dispositivo stesso. L'allarme sul touch-screen deve essere confermato prima di poter continuare con il funzionamento.

Se viene emesso un allarme, sul display touch-screen appare una barra rossa in alto sullo schermo attuale. Sull'area di informazione e di stato operativo (⑦ nella Figura 26 e ⑧ nella Figura 27) appare una campana d'allarme rossa racchiusa da due linee di squillo. Inoltre, un segnale d'allarme acustico emette un suono continuo.

Sotto la barra rossa d'allarme un messaggio a testo scorrevole visualizza un riassunto dell'attuale allarme. Sul lato destro appare un pulsante di snooze che permette di tacitare temporaneamente l'allarme. Se la condizione d'allarme non viene eliminata entro il periodo di snooze, il segnale acustico d'allarme riprende a suonare. La durata della tacitazione temporanea (snooze) può essere scelta nelle impostazioni, come spiegato alla sezione «Allarmi e avvertimenti» a pagina 70.

Toccando il simbolo della campana nell'area di informazione e di stato operativo (2) nella Figura 26 e (3) nella Figura 27) si apre una schermata che presenta tutti i dettagli dell'allarme attualmente attivo, come illustrato nella Figura 37. Si ha la possibilità di scorrere l'elenco e di toccare qualsiasi voce dell'elenco per estendere il testo e visualizzare maggiori dettagli.



Figura 37: Lista degli allarmi

Si ha la possibilità di selezionare l'allarme attivo, spuntando la casella accanto alla voce di allarme. Toccando il pulsante **Conferma** il sistema cerca di cancellare l'allarme dalla lista se la causa dell'allarme è stata eliminata. Dopo la cancellazione di tutti gli allarmi l'icona dello stato ritorna al simbolo del cuore verde.

<u>Errore</u>

Al verificarsi di un malfunzionamento, l'agitatore visualizza un messaggio d'errore e si ferma subito per evitare danni ai campioni e/o al dispositivo stesso. Lo schermo diventa completamente rosso e non sono possibili ulteriori interazioni. Viene visualizzato un messaggio di errore insieme a un codice d'errore, come illustrato nell'esempio riportato nella Figura 38.



Figura 38: Schermata di errore

Procedere come segue per porre rimedio al problema e per cercare di rimettere in funzione il sistema:

- 1. Prendere nota del codice d'errore indicato nella schermata di errore.
- 2. Toccare **Conferma** per tacitare il segnale acustico di allarme.
- 3. Spegnere e riaccendere l'agitatore per riavviarlo.
- Se il messaggio di errore continua a essere presente, contattare il servizio di assistenza ai clienti di Thermo Fisher Scientific e comunicare il codice d'errore indicato sulla schermata di errore.

3. 2. 3. Impostazioni

Il secondo simbolo nella barra di navigazione è il simbolo delle impostazioni. Toccando **Impostazioni** appare la schermata illustrata in basso. La schermata Impostazioni offre un numero di pulsanti superiore a quelli visualizzabili sul display. Per vedere gli altri pulsanti scorrere verso il basso, come illustrato sul lato destro della Figura 39.

<	Shake Lab 1 11:22:24AM Settings		<	Shake Lab 1 11:22:24AM Settings	
۵.	Alarms and Alerts	î	\$₀ (Controls	ÎŶ
Ë	Access Control		邕	🖵 Display	
Ē	Calibration		Ē	Files and Info	
Ļ	Controls	~	Ļ	🗙 Service 🏔	- ~
		()			()

Figura 39: Schermata Impostazioni

Allarmi e avvertimenti

La schermata Allarmi e Avvertimenti permette di determinare come e quando devono essere emessi allarmi e avvertimenti.



Figura 40: Impostazioni per gli allarmi e gli avvertimenti

Volume: Questo regolatore permette di modificare il volume del segnale acustico emesso in caso di allarme o avvertimento. Trascinare il cursore verso sinistra per ridurre il volume o verso destra per aumentarlo, quindi toccare **Salva**. L'impostazione di questo volume riguarda sia gli allarmi sia gli avvertimenti.



Figura 41: Impostazione del volume del segnale acustico per gli allarmi e gli avvertimenti

Allarme di temperatura alta / allarme di temperatura bassa (solo per agitatori a termoregolazione): Queste due opzioni permettono di impostare un valore limite per l'allarme di temperatura alta e bassa rispetto al valore nominale della temperatura (vedere «Impostazione della temperatura» a pagina 65). Appena la temperatura all'interno della camera dell'agitatore scende sotto il valore limite inferiore o supera quello superiore l'agitatore emette un allarme. Scegliere un valore limite per la temperatura e toccare **Salva**.



Figura 42: Impostazione dell'allarme di temperatura alta e bassa

AVVISO Se viene modificato il valore nominale della temperatura, i valori limite per l'allarme si spostano rispettivamente. Spegnimento per temperatura alta / spegnimento per temperatura bassa (solo per agitatori a termoregolazione): Queste due opzioni permettono di impostare un valore limite per lo spegnimento di sicurezza per temperatura alta e bassa rispetto all'attuale valore nominale della temperatura (vedere «Impostazione della temperatura» a pagina 65). Appena la temperatura all'interno della camera dell'agitatore scende sotto il valore limite di temperatura bassa o supera il valore limite di temperatura alta, l'agitatore si spegne automaticamente per proteggere i campioni che si trovano nella camera. Scegliere un valore limite e toccare **Salva**.

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

Figura 43: Impostazione dei valori limite superiore e inferiore per lo spegnimento

AVVISO Se viene modificato il valore nominale della temperatura, i valori limite di errore si spostano rispettivamente.

Durata della tacitazione temporanea (Snooze): La durata della tacitazione temporanea determina per quanto tempo l'allarme viene tacitato dopo aver toccato il pulsante di snooze sulla barra rossa che appare in alto nella finestra principale (vedere «Allarme» a pagina 68). La durata della tacitazione può essere impostata a 5 minuti, 10 minuti o 15 minuti. L'opzione di default è 10 minuti. Toccare il pulsante **Salva** per salvare le modifiche apportate.



Figura 44: Impostazione della durata della tacitazione per gli allarmi

Disattivazione degli avvisi di calibrazione (solo per agitatori a termoregolazione): Questa opzione permette di disattivare gli avvisi che ricordano l'utente a intervalli regolari di calibrare la misurazione della temperatura sull'agitatore (per maggiori informazioni consultare «Calibrazione di temperatura» a pagina 127). Si ha la possibilità di scegliere tra la disattivazione completa degli avvisi di calibrazione o la disattivazione durante l'esecuzione di un programma.

Disable Calibration Notifications	Disable Calibration Notifications while a program is actively running
Figura 45: Disattivazione d	egli avvisi di calibrazione

Controllo dell'accesso

La schermata Controllo all'accesso permette di assegnare un codice di accesso da amministratore per gestire i processi con diritti di accesso limitati.

Per default, l'agitatore si trova in modalità «Senza controllo», cioè non è richiesto l'inserimento di un codice di accesso per accedere e utilizzare il dispositivo. La commutazione sulla modalità «Con controllo dell'accesso» richiede l'inserimento del codice da amministratore per modificare le impostazioni dell'agitatore. Viene visualizzato un pulsante di login nell'angolo in alto a destra dello schermo touch-screen.

Il seguente confronto delle modalità «Senza controllo» e «Con controllo» indica le situazioni in cui è richiesto il login con codice di accesso.

Azione	Codice di accesso richiesto in modalità «Senza controllo»	Codice di accesso richiesto in modalità «Con controllo»
Impostazione dei valori nominali per velocità, durata e temperatura	No	No
Avvio e arresto dell'agitatore	No	No
Calcolo delle orbite	No	No
Confermare allarmi e avvertimenti	No	No
Visualizzare la schermata File & Info	No	No
Visualizzazione dello stato operativo	No	No
Visualizzazione ed esportazione del registro degli eventi e dei diagrammi	No	No
Tacitare temporaneamente gli allarmi	No	No
Eseguire programmi	No	No
Creare, modificare ed eliminare programmi	No	Sì
Modificare le impostazioni del display	No	Sì
Modificare i valori impostati	No	Sì
Modifica delle impostazioni di allarme e avvertimento	No	Sì
Modifica delle impostazioni per il controllo all'accesso	No	Sì
Importazione o esportazione dei programmi	No	Sì
Ripristino delle impostazioni di fabbrica	No	Sì
Effettuazione di una calibrazione	No	Sì
Installazione di un aggiornamento firmware	Sì	Sì

Tabella 52: Necessità di effettuare il login con codice di accesso in modalità «Senza controllo» e «Con controllo»
La Figura 46 illustra la schermata di Controllo dell'accesso in modalità «senza controllo» e «con controllo».



Figura 46: Controllo dell'accesso: Modalità «senza controllo» (a sinistra) e «con controllo» (a destra)

Selezionando **Con controllo** e toccando quindi il pulsante **Salva** il sistema chiede l'inserimento del codice da amministratore prima che la modifica possa essere confermata. Tutte le unità sono consegnate con lo stesso codice di accesso per amministratori preimpostato in fabbrica e riportato nel manuale.

Allo stesso modo, a qualsiasi utente che prova di modificare un'impostazione protetta da codice di accesso viene chiesto l'inserimento del codice da amministratore.

AVVISO Il codice di amministratore preimpostato è «00000».

Il pulsante **Modifica codice di accesso Admin** permette di modificare il codice di accesso preimpostato. Toccando il pulsante **Modifica codice di accesso Admin** il sistema chiede all'utente di inserire prima l'attuale codice di accesso, e poi un'altra volta il nuovo codice di accesso per confermarlo. Il messaggio **Il nuovo codice di accesso**.

Calibrazione

La schermata Calibrazione (solo per agitatori a termoregolazione) apre una serie di schermate che permettono di calibrare l'agitatore in modo da ottenere una regolazione precisa della temperatura. La procedura è descritta in «Calibrazione di temperatura» a pagina 127.

Elementi di comando

La schermata Elementi di comando offre opzioni per la preimpostazione dei parametri operativi di default attivi all'avvio del dispositivo:

Valore nominale di velocità: Il valore nominale di velocità può essere preimpostato a un valore tra 15 e 525 giri/min. Ruotare i quadranti della rotella di selezione per preimpostare la velocità e toccare **Salva**. Per informazioni dettagliate sull'uso delle schermate della velocità nominale consultare la sezione «Impostazione della velocità» a pagina 62 al capitolo 3. 2. 1.

Modalità tempo: L'agitatore può essere usato in modalità continua o temporizzata. In modalità continua l'agitatore può essere fermato manualmente a propria discrezione. In modalità temporizzata l'azionamento dell'agitatore si ferma automaticamente dopo che è trascorso il tempo impostato. Per la modalità temporizzata si ha la possibilità di preimpostare una durata di default e scegliere se durante il funzionamento dell'agitatore l'interfaccia utente deve visualizzare il tempo trascorso oppure il tempo rimanente. Per informazioni dettagliate sull'uso della schermata della modalità temporizzata consultare la sezione «Impostazione della durata del ciclo» a pagina 64 al capitolo 3. 2. 1.

Valore nominale della temperatura (solo per agitatori a termoregolazione): Questa opzione permette di impostare una temperatura di default visualizzata dall'agitatore al momento dell'avvio. Per informazioni dettagliate sull'uso della schermata della temperatura nominale consultare la sezione «Impostazione della temperatura» a pagina 65 al capitolo 3. 2. 1.

Riavvio automatico: Questa funzione riavvia il dispositivo dopo un'interruzione dell'alimentazione che si è verificata durante un normale ciclo di avvio/arresto, durante l'esecuzione di un programma o durante un ciclo di calibrazione della temperatura. Se il Riavvio automatico è impostato su **No**, il dispositivo non sarà riavviato dopo un'interruzione dell'alimentazione.

Display

Le impostazioni del display permettono di modificare diverse opzioni di visualizzazione.

Luminosità: Per regolare il livello di luminosità del display usare il comando a cursore oppure i pulsanti +/-.



Figura 47: Regolazione della luminosità dello schermo

Lingua: Per modificare la lingua di visualizzazione toccare il pulsante Lingua. Selezionare la lingua desiderata ruotando la rotella di selezione e toccare Salva.



Figura 48: Selezione della lingua di visualizzazione

Unità di misura (solo per agitatori a termoregolazione): Toccare Unità di misura per commutare tutti i valori di temperatura visualizzati sulle schermate dell'interfaccia utente tra gradi centigradi °C e gradi Fahrenheit °F.

Units of Measure	°C 🔨
	۴F

Figura 49: Selezione dell'unità di misura per la temperatura

Data: Per impostare la data, toccare il pulsante Data. Ruotare i quadranti per il mese, il giorno e l'anno della rotella di selezione e toccare **Salva**.



Figura 50: Impostazione di mese, data e ora

Formato data: Per impostare il formato data, toccare il pulsante **Formato data**. Toccare il pulsante di opzione per il formato di data desiderato (esempio: con MM/ GG/AAAA la data appare nel formato Aprile 15, 2015) e toccare **Salva**.



Figura 51: Impostazione del formato data

Tempo: Per impostare l'ora e il formato dell'ora, toccare il pulsante **Ora**. Toccare **12ore** o **24ore**, quindi ruotare il quadrante delle ore, dei minuti e AM/FM (solo per il formato a 12 ore) della rotella di selezione e toccare **Salva**.



Figura 52: Impostazione dell'ora e del formato 12 ore o 24 ore

Modo veglia: Toccare **Modalità Veglia** per portare il display dell'agitatore in modalità veglia dopo un periodo di inattività di 15 minuti. In modalità veglia lo schermo visualizza l'invito a **Toccare per attivare**.



Figura 53: Attivazione della modalità veglia

Regione: Per l'impostazione della regione in cui il dispositivo viene usato toccare il pulsante **Regione**. Toccare il campo di testo Città, Paese e iniziare a digitare i primi caratteri del nome della città. Dopo l'inserimento dei primi tre caratteri l'interfaccia utente inizia a suggerire possibili corrispondenze. Accettare una corrispondenza o continuare a digitare il nome intero, poi toccare **Salva** sulla tastiera su schermo.

< *	Sh 11: Re	ake La 22:24/ gion	b 1 AM			0					< ♠	Shake Lab 1 11:22:24AM Region	\bigcirc	
≎ 。 ₿	Cit	y, Cour	ntry:	Enter the	city and i	country (closest to	your tin	te zone		¢, (≝	City, Country:	Enter the city and country closest to your time zone New York, United States	
Q	w	E	R	т	Y	U	1	0	Ρ	€	Ē			
A	S	D	F	G	Н	J	к	L	Sa	ve				
1	Z	x	С	V	В	N	м			1				
?!1:	23								-	W			Save	()

Figura 54: Imposta Regione

Nome dell'unità: Per inserire o modificare il nome dell'agitatore, toccare il pulsante Nome dell'unità. Toccare il campo di testo Nome dell'unità e iniziare a digitare il nome. Alla fine toccare il pulsante **Salva** sulla tastiera su schermo.

< ♠	Sh 11 Un	ake La :22:24/ it Nam	b 1 AM ie			2					< *	Shake Lab 1 11:22:24AM Unit Name	\bigcirc	
٥.	U	nit Na	me	Sha	ake La	ab 1					¢.			
Q	W	E	R	Т	Y	U	1	0	Ρ			Unit Name:	Shake Lab 1	
А	S	D	F	G	Н	J	K	L	Sa	ave				
1	Z	X	С	V	В	Ν	М	0		1			Save	0
?!1:	23							-	-	÷.				U

Figura 55: Assegnazione di un nome all'agitatore

Personalizzazione del menu: Toccare il pulsante **Personalizza menu** per personalizzare due simboli in fondo alla barra di navigazione principale. Trascinare il simbolo desiderato dall'area della schermata principale sul simbolo che si desidera sostituire. Toccare **Salva** per confermare.

AVVISO Il pulsante **Resetta menu** permette di ripristinare in qualsiasi momento le impostazioni di fabbrica per la barra di navigazione.



Figura 56: Personalizzazione della barra di navigazione

File & Info

Questa schermata visualizza il numero di serie, il software applicativo e la versione dell'interfaccia utente, il firmware e la versione del controllore principale dell'agitatore, il file parametri e la versione nonché l'attuale sistema operativo.

Il pulsante **Reset ai valori di fabbrica** permette inoltre di ripristinare le impostazioni di fabbrica per l'agitatore. Il ripristino delle impostazioni di fabbrica richiede l'inserimento del codice da amministratore e cancella tutte le impostazioni effettuate attraverso l'interfaccia utente grafica. Il registro degli eventi non viene cancellato dal ripristino delle impostazioni di fabbrica.

< ☆ ≝	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	 < Sha 11:2 Files ♣< < 	ke Lab 1 12:24AM a 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	>	
.	Factory Reset	.	Firmware	>	Ŷ
	0				()

Figura 57: File & Info

Se si scorre la schermata verso il basso appare il pulsante **Firmware** illustrato sul lato destro della Figura 57. Questo pulsante può essere toccato per installare un nuovo firmware per l'agitatore. Per le istruzioni dettagliate sull'installazione del nuovo firmware consultare «4. 7. Installazione del firmware» a pagina 131.

AVVISO Un nuovo firmware deve essere fornito da un tecnico di assistenza post vendita autorizzato.

Manutenzione

L'accesso alle impostazioni di servizio è limitato ai tecnici di assistenza post vendita autorizzati.

3. 2. 4. Programmi

Toccare il simbolo **Programmi** nella barra di navigazione per visualizzare la lista dei programmo o per creare un nuovo programma. Si ha la possibilità di creare, modificare, eliminare, importare ed esportare un programma. Nella Figura 58 è illustrata la schermata di un dispositivo nuovo (a sinistra) e di un dispositivo già utilizzato da tempo (a destra). Il dispositivo già in uso visualizza una lista di programmi creati dall'utente.



Figura 58: Programmi

Ognuna delle voci nella lista di programmi dispone di un simbolo di occhio () che apre una schermata di vista rapida con dettagli sul programma e un simbolo di penna / che apre la schermata di editing per il programma.

Impostazione delle preferenze per tutti i programmi

Dalla schermata Programmi, con il pulsante **Impost. avanzate** si accede alla schermata delle **Impostazioni avanzate**. Questa schermata permette di determinare cosa succede all'avvio di un programma. Queste preferenze sono applicate a tutti i programmi che si creano.



Figura 59: Impostazione delle preferenze per i programmi

1. Toccare il pulsante Impost. avanzate.

- 2. Scegliere un'opzione per **Mostra timer** per determinare in che modo viene visualizzato il temporizzatore del programma:
 - » Tempo trascorso: tempo trascorso, cioè per quanto tempo l'agitatore sta già funzionando dal momento in cui è stato premuto il pulsante di avvio, oppure
 - » Tempo rimanente: tempo residuo, cioè per quanto tempo l'agitatore continuerà a funzionare fino al trascorrere del tempo impostato.
- Se si desidera che l'agitatore parte subito dopo la scelta di un programma, attivare l'opzione Avvia agitatore al caricamento programma. L'impostazione predefinita prevede che si deve toccare il pulsante Avvio per eseguire il programma.

AVVISO L'opzione **Avvia agitatore al caricamento programma** funziona solo con la cappa dell'agitatore chiusa. Appena viene chiusa la cappa, il ciclo parte automaticamente.

4. Toccare il pulsante **Salva**.

Creazione di un programma

Possono essere creati e salvati un totale di 99 programmi.

- 1. Toccare il pulsante Crea Nuovo.
- 2. Inserire il nome del programma.



Figura 60: Creare un programma

3. Inserire la temperatura (solo per agitatori a termoregolazione), la velocità e la durata del programma in ore e minuti.

Per disattivare la regolazione della temperatura su un agitatore a termoregolazione lasciare vuoto il campo**Temp**.

- 4. Per aggiungere uno step supplementare al programma, scorrere giù e toccare il pulsante **Aggiungi step**.
- 5. Toccare il pulsante **Salva** per salvare il programma.

Modifica di un programma

- 1. Toccare il simbolo della penna 🖍 accanto al programma che si desidera modificare.
- 2. Modificare i campi desiderati. Toccare il pulsante **Salva**. Il programma viene salvato con le modifiche apportate.
- Si ha la possibilità di aggiungere uno step, selezionando Aggiungi step in fondo alla schermata. Scorrere la schermata per vedere il pulsante Aggiungi step, se sono presenti più di 3 step.

AVVISO Se durante una procedura di agitazione composta di diversi step si verifica un messaggio d'errore o d'allarme, l'agitatore si ferma automaticamente. Se viene visualizzato un messaggio d'avvertimento, la procedura di agitazione continua.



Figura 61: Aggiungere un passo

Eliminazione di un programma

- 1. Toccare il simbolo della penna accanto al programma che si desidera modificare.
- 2. Toccare il pulsante **Elimina**. Appare una finestra che chiede una conferma.



Figura 62: Eliminare un programma

3. Toccare il pulsante **Elimina**. Il programma selezionato viene eliminato.

Esecuzione di un programma

- 1. Per eseguire un programma esistente, selezionare il programma che si desidera eseguire.
- Per accedere a una vista rapida del programma, toccare il simbolo dell'occhio
 accanto al programma.
- 3. Appare una finestra con una visione sintetica del programma.



Figura 63: Vista rapida del programma

- Toccare il pulsante Ciclo per caricare il programma sulla schermata Home.
 In base alle impostazioni di Avvio automatico scelte, il programma viene avviato subito oppure con l'azionamento del pulsante Avvio sulla schermata principale.
- 5. Dopo il completamento del programma appare la seguente finestra. Toccare il pulsante **OK**.



Figura 64: Programma completato

Importa programmi

I programmi creati su un agitatore possono essere importati su un altro agitatore. A tale scopo il dispositivo USB deve essere inserito nella porta USB dell'agitatore.

La schermata seguente visualizza tutti i programmi identificati sul dispositivo USB.



Figura 65: Selezionare i programmi da importare

Selezionare i programmi che si desiderano importare. Toccare il pulsante **Importazione**.

Dopo l'importazione riuscita dei programmi appare la finestra riportata di seguito. Toccare il pulsante **OK**. Adesso la chiavetta USB può essere rimossa.



Figura 66: Importazione programma completata

Esportazione dei programmi

I programmi possono essere esportati da un agitatore ad un altro agitatore. Assicurarsi che sia collegata una chiavetta USB.

Se è collegata una chiavetta USB, viene visualizzata la seguente schermata:



Figura 67: Selezionare i programmi da esportare

Selezionare i programmi che si desiderano esportare. Toccare **Esporta**.

Al completamento dell'esportazione viene visualizzata la seguente finestra. Toccare il pulsante **OK**. Adesso la chiavetta USB può essere rimossa.



Figura 68: Esportazione programma completata

Registro degli eventi

Il terzo pulsante nella barra di navigazione è il registro degli eventi che contiene un record degli eventi utente e sistema. La schermata Registro eventi viene visualizzata toccando il simbolo Registro Eventi nella barra dei menu, come illustrato nella Figura 69.



Figura 69: Registro degli eventi

Questa schermata visualizza una lista degli eventi recenti, insieme alla data e all'ora per ogni evento.

Selezionando una voce dalla lista è possibile visualizzare informazioni aggiuntive per il singolo evento.

Gli eventi possono essere filtrati secondo le seguenti categorie:

- allarmi
- avvertimenti
- modifiche apportate alle impostazioni
- azioni di avvio/arresto
- cicli di programma
- azioni di apertura cappa (solo per agitatori a termoregolazione)

Se viene selezionato un filtro, il pulsante Visualizza tutto a destra cambia in Filtro ON, come illustrato nella Figura 70.

	Shake Lab 1 11:22:24AM Event Log	\checkmark)		
A	Date/Time	Event	Filter ON	\mathbf{v}	^
Q .	5/12/15 11:34:06AM	Setpoint Change		$\mathbf{\sim}$	
<u>é</u>	5/12/15 11:34:06AM	High Temp Alarm		\checkmark	
	5/12/15 11:34:06AM	Unit Stopped		\checkmark	
Ļ	5/12/15 11-24-06AM	Unit Started		\checkmark	~
	Export	Log	View Cha	rt	Ó

Figura 70: Registro degli eventi con filtro

Esportazione del registro degli eventi

1. Selezionare gli eventi da esportare dalla lista a discesa **Eventi da esportare**. Selezionare il formato di esportazione del registro o report.



Figura 71: Selezionare un registro degli eventi da esportare

2. Può essere selezionato un periodo predefinito o inserito un campo data personalizzato.

AVVISO Possono essere esportati solo eventi verificatesi negli ultimi sei mesi.

 L'esportazione può essere effettuata con l'ausilio di una chiavetta USB. Deve essere inserita una chiavetta USB per salvare il registro o il report. Toccare il pulsante **Esporta registro** per scaricare il registro o il report.



Figura 72: Inserire la chiavetta USB per l'esportazione

<u>Diagrammi</u>

I diagrammi visualizzano grafici con i dati di velocità o temperatura (solo per agitatori a termoregolazione) nel tempo. L'asse X riporta il tempo e l'asse Y visualizza la velocità o la temperatura.

Il diagramma di temperatura permette di registrare i valori per la temperatura nominale, effettiva e per la temperatura ambiente nel tempo. Il pulsante **Modifica** permette di scegliere quale delle tre curve si vuole includere nel diagramma della temperatura. In più, è possibile specificare un campo di visualizzazione di 1 giorno, 7 giorni o di una serie di giorni specificati dall'utente, per le intere 24 ore oppure per le ore del giorno selezionate.

Il pulsante **Aggiorna** aggiorna la visualizzazione del diagramma sulla base delle impostazioni effettuate e visualizza i nuovi dati registrati dall'ultima volta che il diagramma è stato caricato o aggiornato.

AVVISO In caso di un allarme temperatura, la temperatura visualizzata sull'interfaccia grafica utente passa immediatamente alla temperatura effettiva dell'aria all'interno della camera dell'agitatore per proteggere al meglio i campioni. Ciò potrebbe comportare un salto della temperatura della camera visualizzata nella funzione grafico.



Figura 73: Diagrammi di velocità e temperatura

Si ha la possibilità di inserire un dispositivo USB nella porta USB dell'agitatore e toccare **Esporta dati diagramma** per scaricare i dati del diagramma di velocità e temperatura.

Modifica del diagramma

1. Toccare **Modifica** per avviare la modifica del diagramma.



Figura 74: Modifica del diagramma: Selezione della temperatura o della velocità da registrare

2. Toccare **Temp** o **Velocità** per scegliere il diagramma da visualizzare.

- 3. Scorrere giù nella schermata per visualizzare le opzioni Campo Data.
- 4. Toccare 1 giorno, 7 giorni o **Personalizza** per specificare un campo data personalizzato.
- 5. Ruotare i quadranti della rotella di selezione per scegliere la data di inizio e toccare il pulsante **Imposta Data iniziale**.
- Ruotare i quadranti della rotella di selezione per scegliere la data di fine e toccare il pulsante **Imposta Data finale**.



Figura 75: Selezione del campo data per il diagramma

- 7. Scorrere giù nella schermata per visualizzare le opzioni di Campo Ora.
- 8. Per effettuare la selezione toccare 24ore al giorno oppure **Personalizza** per registrare un periodo di tempo specifico e
 - a. Ruotare i quadranti della rotella di selezione per scegliere l'ora iniziale quotidiana e toccare il pulsante **Imposta Ora iniziale**.
 - b. Ruotare i quadranti della rotella di selezione per scegliere l'ora finale quotidiana e toccare il pulsante **Imposta Ora finale**.

AVVISO Con la scelta di un'ora iniziale e un'ora finale si decide di raffigurare i dati per un periodo esatto tra due punti temporali. Se si lavora con l'opzione 24 ore al giorno, la registrazione inizia alle ore 00:00 e termina alle ore 23:59 di ogni giorno.



Figura 76: Selezione delle ore del giorno per la registrazione

9. Toccare il pulsante **Salva** per salvare le modifiche apportate o le impostazioni personalizzate.

3.3. Accessori



Lesioni e rischio biologico causato da recipienti rotti. Accessori non installati correttamente possono comportare la rottura di vetro e la fuoriuscita dei campioni.

ATTENZIONE

Assicurarsi che gli accessori siano installati accuratamente con l'ausilio degli attrezzi e delle viti corretti.

Assicurarsi che gli accessori siano adatti alla piattaforma. Utilizzare i recipienti sempre con gli accessori adatti alla rispettiva dimensione.



ATTENZIONE

Ferite da taglio causate da spigoli taglienti.

Fare attenzione durante la manipolazione delle piattaforme e di altri accessori.



Possibile schiacciamento delle dita causato da una piattaforma in movimento.

ATTENZIONE

Non mettere mai le dita su o sotto una piattaforma in movimento.

AVVISO L'operatore stesso è responsabile della corretta installazione degli accessori.

Le procedure d'installazione descritte si applicano a tutti i modelli di agitatore elencati nelle presente istruzioni, fino a diversa specificazione.

Utilizzare sempre gli attrezzi corretti, in particolare gli attrezzi forniti insieme all'accessorio. Qualora si dovessero perdere alcuni degli articoli, i kit di ricambi elencati potranno essere riordinati («Accessori» a pagina 18). Non utilizzare attrezzi diversi da quelli specificati o viti diverse da quelle fornite.

3. 3. 1. Installazione della piattaforma

ATTENZIONE Una lista completa delle piattaforme compatibili per ogni agitatore si trova alla sezione «1. 2. 1. Piattaforme» a pagina 19. I modelli di agitatori a termoregolazione devono essere utilizzati solo con le piattaforme preinstallate in fabbrica.

ATTENZIONE Ferite da taglio causate da spigoli taglienti. Per smontare una piattaforma, mettere le mani sotto la piattaforma. Non sollevare la piattaforma afferrando i supporti montati sulla piattaforma. Fare attenzione durante la manipolazione delle piattaforme e di altri accessori.

ATTENZIONE Nell'agitatore Solaris 4000 I / 4000 R sussiste il rischio di schiacciamento della mano o delle dita con la piattaforma 11x14 che si muove. Non usare la piattaforma 11x14 nell'agitatore Solaris 4000 I / 4000 R.

ATTENZIONE Utilizzare solo le viti e gli attrezzi di serraggio forniti insieme alla piattaforma. L'utilizzo di viti diverse o di attrezzi di serraggio sbagliati comporta un'installazione non corretta e possibili danni all'agitatore e agli accessori.

Installare una piattaforma sempre con tutte le viti.

Utilizzare sempre l'attrezzo di serraggio idoneo:

- Attrezzo di fissaggio da 3/16" (GT530066) per tutte le piattaforme dell'agitatore Solaris 2000 e Solaris 2000 I/R e 4000 I/R
- Attrezzo di fissaggio da 7/32" (GT530080) per tutte le piattaforme dell'agitatore Solaris 4000

Piattaforme universali

Ogni agitatore della serie Solaris 2000 e 4000 viene fornito con una piattaforma universale, viti e un attrezzo di fissaggio. Si ha la possibilità di acquistare ulteriori piattaforme per la specifica applicazione. Una lista completa delle piattaforme compatibili per ogni agitatore si trova alla sezione «1. 2. 1. Piattaforme» a pagina 19. Gli agitatori 2000 I, 2000 R, 4000 I e 4000 R hanno la piattaforma già preinstallata nell'agitatore. Né la piattaforma né le viti sono disponibili come componenti separati.



Figura 78: Piattaforma universale installata su un agitatore Solaris 2000 (a sinistra) e Solaris 4000 (a destra)

1. Collocare la piattaforma con cautela in posizione orizzontale sopra l'agitatore e sui relativi punti di montaggio.

Le piattaforme per gli agitatori Solaris 2000 hanno 3 fori di montaggio. Controllare i campioni, le condizioni ambientali e/o modificare le impostazioni. Platforms for Solaris 4000 shaker have 4 mounting holes.

Le piattaforme per gli agitatori Solaris 2000 l/R e 4000 l/R hanno 3 fori di montaggio.
 Serrare le viti delicatamente per fissare la piattaforma all'agitatore. Appena l'attrezzo di serraggio inizia a curvarsi, interrompere il serraggio della vite.

Piattaforme a due livelli

AVVISO Piattaforme a due livelli non sono compatibili con l'agitatore Solaris 2000 I/R e 4000 I/R.



Figura 79: Piattaforma a due livelli

Collegare la piattaforma superiore e quella inferiore, montando i 4 montanti, uno ad ogni angolo. Fissare i montanti con le corrispondenti viti alla piattaforma superiore e alla piattaforma inferiore.

Serrare le viti delicatamente per collegare i montanti e le piattaforme. Appena l'attrezzo di serraggio inizia a curvarsi, interrompere il serraggio della vite.

La piattaforma inferiore è quella dotata dei fori di montaggio per il collegamento all'agitatore. Durante l'assemblaggio assicurarsi di collocare la piattaforma inferiore in basso.

Accessori

AVVISO Prima di posizionare dei recipienti sul gruppo piattaforma, effettuare un ultimo controllo per assicurarsi che il gruppo piattaforma non oscilli.



Figura 80: Piattaforma a due livelli su un agitatore Solaris 4000

1. Collocare la piattaforma a due livelli assemblata con cautela in posizione orizzontale sull'agitatore e sulla piastra di montaggio piattaforma.

Le piattaforme per Solaris 2000 hanno 3 fori di montaggio, le piattaforme per Solaris 4000 ne hanno 4.

2. Serrare le viti delicatamente per fissare la piattaforma all'agitatore. Appena l'attrezzo di serraggio inizia a curvarsi, interrompere il serraggio della vite.

3. 3. 2. Installazione dei supporti porta matracci e dei recipienti

ATTENZIONE Rischio biologico dovuto a recipienti rotti o che perdono. Accessori non installati correttamente possono comportare la fuoriuscita dei campioni. Assicurarsi che gli accessori siano installati accuratamente con l'ausilio degli attrezzi e delle viti corretti. L'installazione richiede un cacciavite (PH2) con una lunghezza di stelo di 150 mm / 6 pollici. (Numero articolo 75004131). Assicurarsi che gli accessori siano adatti alla piattaforma. Utilizzare i recipienti sempre con gli accessori adatti alla rispettiva dimensione. I recipienti devono essere intatti e installati accuratamente.

ATTENZIONE Ferite da taglio causate da spigoli taglienti. Fare attenzione durante la manipolazione delle piattaforme e di altri accessori.

ATTENZIONE Un fissaggio sicuro può essere garantito solo in combinazione con le viti fornite. Inserire le viti fornite in tutti i fori di montaggio previsti per il supporto.

AVVISO Se le viti sono troppo lunghe ciò può pregiudicare la stabilità e la funzionalità della piattaforma intercambiabile. Fare attenzione a utilizzare per il fissaggio sempre solo le viti comprese nella fornitura.

Supporti porta beute

Ogni supporto porta beute è composto di una clip porta beute/pallone, una o due molle (in base alle dimensioni del supporto porta beute/flacone) e viti per il montaggio alla piattaforma. Utilizzare solo le viti fornite insieme al supporto porta beute.





① Bussola a molla ② Molla ③ Viti ④ Clip ⑤ Molla

Figura 81: Supporto porta beute con 2 molle

- 1. Se necessario, fissare la molla alle clip come illustrato nella figura.
- Le bussole a molla in gomma sono posizionate tra le clip, come illustrato nella figura. Alcuni supporti porta beute usano due molle. La seconda molla viene installata intorno alla base del gruppo clip dopo l'avvenuto montaggio sulla piattaforma.
- 3. Con l'ausilio delle viti in dotazione, fissare il gruppo clip sulla piattaforma.

Recipiente

 Posizionare il recipiente desiderato con cautela nel supporto porta beute, tirando la molla quanto basta in maniera sufficiente per da consentire l'introduzione della base del recipiente nel supporto porta beute. Abbassare il recipiente delicatamente fino alla posizione corretta e bloccarlo nel fondo più largo del supporto porta beute. La molla si stringe sul collo del recipiente e lo mantiene in posizione sicura.



2. Assicurarsi che tutti i recipienti siano fissati accuratamente prima di accendere il dispositivo.

Ove possibile, il recipiente dovrebbe essere chiuso con un tappo per impedire che le sostanze possano essere catapultate fuori durante la miscelazione.

3. 3. 3. Installazione dei supporti quadrati

ATTENZIONE Rischio biologico dovuto a recipienti rotti o che perdono. Accessori non installati correttamente possono comportare la fuoriuscita dei campioni. Assicurarsi che gli accessori siano installati accuratamente con l'ausilio degli attrezzi e delle viti corretti. L'installazione richiede un cacciavite (PH2) con una lunghezza di stelo di 150 mm / 6 pollici. (Numero articolo 75004131). Assicurarsi che gli accessori siano adatti alla piattaforma. Utilizzare i recipienti sempre con gli accessori adatti alla rispettiva dimensione. I recipienti devono essere intatti e installati accuratamente.

ATTENZIONE Ferite da taglio causate da spigoli taglienti. Fare attenzione durante la manipolazione delle piattaforme e di altri accessori.

Dettagli del supporto quadrato

Ogni supporto quadrato è composto di una piastra di fondo con fori di montaggio e quattro pareti collegate a molla. Inoltre, sono comprese le viti per il montaggio del supporto sulla piattaforma. Utilizzare solo le viti fornite insieme al supporto porta beute.



① Fori di montaggio ② Molla ③ Parete collegata a molla ④ Viti

Figura 82: Dettagli del supporto quadrato

- 1. Se necessario, agganciare le molle 2 nelle pareti 3, come illustrato nella Figura 82.
- 2. Posizionare il supporto assemblato sulla piattaforma e allineare i fori di montaggio ①.
- 3. Fissare il supporto assemblato alla piattaforma con le viti ④ fornite.
- 4. Inserire il recipiente quadrato dall'alto.
- 5. Assicurarsi che il recipiente quadrato poggi sul fondo e sia posizionato saldamente nel supporto.

3. 3. 4. Installazione dei rack portaprovette

ATTENZIONE Ferite da taglio causate da spigoli taglienti. Fare attenzione durante la manipolazione delle piattaforme e di altri accessori.

Ogni rack portaprovette è composto di due staffe di montaggio (in filo metallico o lamiera) con un pomello di fissaggio e due viti zigrinate per staffa che consentono di montare il supporto alla piattaforma. Ogni unità di rack portaprovette è consegnata preassemblata con un rack portaprovette e inserti in materiale espanso. Utilizzare sempre le viti fornite insieme al rack.



Figura 83: Unità di rack portaprovette con staffe di montaggio in filo metallico



- Rack portaprovette
 Inserti in materiale espanso
- 3 Pomello di fissaggio
- ④ Staffa di montaggio

Figura 84: Unità di rack portaprovette con staffe di montaggio in lamiera

- 1. Montare le staffe di montaggio per mezzo delle viti zigrinate sulla piattaforma. Serrare a mano le viti zigrinate.
- 2. Posizionare il rack portaprovette sulle staffe di montaggio.
- Regolare l'angolo richiesto con l'ausilio dei pomelli di fissaggio. Serrare a mano i pomelli di fissaggio.

3. 3. 5. Installazione dei supporti ad angolo regolabile per rack portaprovette

ATTENZIONE Ferite da taglio causate da spigoli taglienti. Fare attenzione durante la manipolazione delle piattaforme e di altri accessori.

I supporti per rack portaprovette e i rack portaprovette sono disponibili in varie dimensioni, come riportato al capitolo «1. 2. 5. Supporti ad angolo regolabile per rack portaprovette» a pagina 29. Tutti i supporti per rack portaprovette sono inclinabili a sette posizioni, ruotando da 0° in entrambe le direzioni e bloccandosi nelle posizioni di 15°, 30° e 45°.

Utilizzare sempre le viti fornite insieme al supporto per rack.



Figura 85: Supporto ad angolo regolabile per rack portaprovette con rack portaprovette installato

- 1. Montare i supporti per rack ① sulla piattaforma dell'agitatore, usando le viti fornite.
- Aprire leggermente le linguette di metallo 3 su entrambi i lati del letto basculante del supporto 6 e inserire il rack portaprovette 2.
- 3. Bloccare il rack portaprovette 2 rilasciando con cautela le linguette 3.
- 4. Tirare i perni di bloccaggio (5) su entrambi i lati del supporto verso l'esterno. I perni vengono sbloccati girando la testa di 1/4 di giro.
- Riposizionare il perno di bloccaggio (5) nel foro di posizionamento (4) e bloccare il perno girando la testa di 1/4 di giro.
- 7. Per rimuovere il rack aprire le linguette di metallo 3 su entrambi i lati del letto basculante del supporto 6 ed estrarre il rack portaprovette 0.

3. 3. 6. Installazione dei supporti per micropiastra / piastra DeepWell

ATTENZIONE Ferite da taglio causate da spigoli taglienti. Fare attenzione durante la manipolazione delle piattaforme e di altri accessori.



 Piastra DeepWell
 Vite
 Supporto per micropiastra

Figura 86: Unità di supporto per micropiastra / piastra DeepWell

- 1. Posizionare il supporto per micropiastra sulla piattaforma.
- Montare il supporto per micropiastra con l'ausilio delle viti fornite insieme al set micropiastra/piastra deepwell alla piattaforma.
- 3. Inserire la micropiastra o piastra deepwell nel supporto.
- 4. Assicurarsi che la micropiastra o la piastra DeepWell sia ferma in posizione, alzandola delicatamente.

3. 3. 7. Installazione dei supporti regolabili per recipienti

ATTENZIONE Ferite da taglio causate da spigoli taglienti. Fare attenzione durante la manipolazione delle piattaforme e di altri accessori.





① Fori di montaggio ② Barre regolabili ③ Pomelli di fissaggio ④ Parete laterale del supporto ⑤ Tappetino antiscivolo ⑥ Viti

Figura 87: Unità di supporto per recipienti a regolazione semplice

Accessori





 ① Fori di montaggio ② Barre regolabili ③ Pomelli di fissaggio ④ Parete laterale del supporto ⑤ Tappetino antiscivolo ⑥ Viti

Figura 88: Unità di supporto per recipienti a doppia regolazione in altezza

- 1. Sollevare il tappetino antiscivolo per esporre i fori di montaggio (posizione 1) nella Figura 88).
- 2. Posizionare il supporto regolabile per recipienti sulla piattaforma e allineare i fori di montaggio.
- 3. Fissare il supporto regolabile per recipienti sulla piattaforma, usando le viti (6) fornite insieme al supporto.
- 4. Posizionare il tappetino antiscivolo (5) sul fondo del supporto regolabile per recipienti.
- 5. Posizionare i recipienti sul tappetino antiscivolo (5).
- 6. Allentare i pomelli di fissaggio 3.
- 7. Regolare le posizioni delle barre regolabili (2) in modo da fissare accuratamente i recipienti.
- 8. Riserrare i pomelli di fissaggio 3.
- 9. Assicurarsi che i recipienti siano bloccati saldamente, cercando di sollevarli leggermente.

3. 3. 8. Installazione del rack porta becher

ATTENZIONE Ferite da taglio causate da spigoli taglienti. Fare attenzione durante la manipolazione delle piattaforme e di altri accessori.



Figura 89: Unità di rack porta becher

- 1. Posizionare la piastra di fondo 2 sulla piattaforma.
- 2. Posizionare il rack porta becher ① sopra la piastra di fondo.
- 3. Allineare i fori di montaggio con la piattaforma.
- 4. Fissare l'unità di rack porta becher con le viti 3 fornite insieme al prodotto.

3. 3. 9. Installazione del supporto per imbuto separatore

ATTENZIONE Ferite da taglio causate da spigoli taglienti. Fare attenzione durante la manipolazione delle piattaforme e di altri accessori.

AVVISO Fare attenzione a utilizzare le viti corrette. Vengono forniti due diversi tipi di viti. Le viti più lunghe con il frenafiletti rosso vanno usati per il fissaggio della staffa verticale. Le viti più corte senza frenafiletti sono utilizzate solo per l'installazione del supporto per imbuto separatore alla piattaforma.

Il supporto per imbuto separatore può essere montato con due orientamenti base alla piattaforma dell'agitatore: in orizzontale e in verticale.



Figura 90: Posizioni di montaggio del supporto per imbuto separatore

La posizione orizzontale ③ nella Figura 90 può essere regolata in altezza per ottenere l'orientamento inclinato illustrato in ④. In più, il supporto per imbuto separatore può essere combinato con la staffa verticale (disponibile come accessorio) per creare i posizionamenti illustrati in ① e ②. Il supporto può essere installato anche con il rubinetto dell'imbuto rivolto verso il basso ① o al contrario per ottenere la configurazione sottosopra illustrata in ②.

Preparazione dell'installazione verticale

La Figura 91 illustra il supporto per imbuto separatore con la staffa verticale (a) accessoria (b) uno accanto all'altro.



Figura 91: Supporto per imbuto separatore (a sinistra) con staffa verticale accessoria (a destra)

- 1. Fissare il supporto per imbuto separatore sul lato lungo della staffa verticale, come illustrato in ① e ② della Figura 90.
- Allineare il supporto per imbuto come descritto di seguito con i quattro fori di montaggio (2) nella Figura 91) della staffa verticale:
 - a. Usare i fori di montaggio per il montaggio verticale standard (3) nella Figura
 91) per un'installazione con il rubinetto rivolto verso il basso.
 - b. Usare i fori di montaggio per l'installazione verticale sottosopra (

 nella
 Figura 91) per l'installazione con il rubinetto rivolto verso l'alto.
 - c. Fissare il supporto per imbuto separatore con le viti fornite (5 nella Figura 91) alla staffa verticale.

Installazione sulla piattaforma

- 1. Posizionare l'unità del supporto per imbuto separatore sulla piattaforma.
- 2. Allineare i fori di montaggio con la piattaforma.
- 3. Fissare il supporto per imbuto separatore alla piattaforma. Usare le viti fornite insieme al supporto per imbuto separatore (6 nella Figura 91 e Figura 92).

Installazione dell'imbuto



Portaimbuto posteriore
Portaimbuto anteriore
Pomello di fissaggio, regolazione in altezza
Pomello di fissaggio, regolazione in lunghezza
Viti senza frenafiletti

1) Imbuto separatore

Figura 92: Unità del supporto per imbuto separatore

- 1. Allentare il pomello di regolazione in lunghezza (5) e scorrere il portaimbuto anteriore (3) sulla massima posizione frontale.
- 2. Inserire l'estremità posteriore dell'imbuto ① nel portaimbuto posteriore ②.
- 3. Spostare il portaimbuto anteriore ③ indietro e posizionare il collo dell'imbuto nell'apertura del portaimbuto anteriore ③.
- Regolare il portaimbuto anteriore ③ a un'altezza conveniente e bloccare la posizione serrando il pomello di regolazione in altezza ④. Con la variazione dell'altezza l'imbuto separatore può essere portato in una posizione inclinata.
- Bloccare il portaimbuto anteriore 3 serrando il pomello di regolazione in lunghezza 6.

3. 3. 10. Installazione del supporto multiuso

ATTENZIONE Ferite da taglio causate da spigoli taglienti. Fare attenzione durante la manipolazione delle piattaforme e di altri accessori.



Figura 93: Unità del supporto multiuso

Per l'installazione procedere come segue:

- 1. Posizionare il supporto multiuso 2 sulla piattaforma.
- 2. Allineare i fori di montaggio con la piattaforma.
- 3. Fissare il supporto multiuso 2 con le viti 3 fornite insieme al supporto multiuso.
- 4. Posizionare il tappetino di gomma ① sul supporto ②.
- 5. Posizionare le bottiglie sul supporto e fissarle con gli elastici O-ring @ forniti, come illustrato nella Figura 94.

AVVISO L'uso degli elastici O-ring è raccomandato per velocità superiori a 100 giri/min.



Figura 94: Supporto multiuso con contenitori fissati e elastici O-Ring

106 / 141

AVVISO Fare attenzione a non tendere eccessivamente gli elastici O-Ring durante il fissaggio dei contenitori. Mentre si tira il singolo elastico durante l'installazione si noterà un punto morto nell'elasticità oltre il quale gli elastici O-Ring non possono più essere tirati. Appena raggiunto questo punto, non continuare a tirare l'elastico per evitare che si strappi.

3. 3. 11. Installazione e uso del tappetino antiscivolo

I tappetini antiscivolo sono idonei per contenitori con una base larga e piatta, come nei matracci di Erlenmeyer e Fernbach, nelle bottiglie recipienti, micropiastre, piastre Petri, fiasche per coltura cellulare, matracci volumetrici e becher. In linea di principio possono essere utilizzati tutti i recipienti che reggono in piedi, ma con la diminuzione del rapporto tra base e altezza diminuisce anche la velocità di agitazione raggiungibile. Contenitori alti con fondo piccolo sono meno adatti, come illustrato nella Figura 95. Matracci di Erlenmeyer grandi (ad es. da 3000 ml) aderiscono meglio rispetto a quelli più piccoli (ad es. da 100 ml).



Figura 95: Tappetino antiscivolo: forme di matracci adatte

Sono disponibili due diversi tipi di tappetino antiscivolo:

- Tappetino antiscivolo a bassa adesione (trasparente) 75004111 e 75004117
 - » gestione facile grazie al basso potere aderente
 - » particolarmente adatto per becher
 - » idoneo per tutti i tipi di recipienti fino a una velocità massima di 200 giri/ min. (applicazioni per la crescita di colture tissutali)
 - Tappetino antiscivolo a alta adesione (nero trasparente) 75004126 e 75004127
 - » fissaggio affidabile dei recipienti grazie all'elevato potere aderente
 - » adatto per tutti i recipienti
 - » velocità massima di 400 giri/min. (fino a Erlenmeyer da 500 ml)

<u>Attrezzi</u>

Il rullo applicatore illustrato nella Figura 96 viene fornito come accessorio per i tappetini antiscivolo. Serve per aiutare a premere il tappeto antiscivolo sulla piattaforma.



Figura 96: Rullo applicatore per tappeto antiscivolo

Posizionamento del tappetino asciutto

- 1. Assicurarsi che la piattaforma dell'agitatore sia pulita e asciutta.
- 2. Rimuovere la pellicola protettiva da un lato del tappetino antiscivolo.
- Incollare il tappetino con la parte adesiva rivolta verso il basso saldamente sulla piattaforma.
- 4. Rimuovere la pellicola protettiva superiore.

Posizionamento del tappetino bagnato

- 1. Rimuovere la pellicola protettiva da entrambi i lati del tappetino antiscivolo.
- 2. Immergere il tappetino in acqua o sciacquarlo accuratamente.
- 3. Fare sgocciolare il tappetino.
- 4. Stendere il tappetino umido sulla piattaforma pulita e posizionarlo come richiesto.
- 5. Fare asciugare il tappetino per 24 ore sulla piattaforma prima di usarlo.

Considerazioni riguardanti le proprietà dei recipienti

La velocità massima dipende dal tipo di recipiente, che si distingue per dimensione, forma, forma del fondo, materiale, rugosità superficiale e pulizia. Questi parametri devono essere valutati per determinare il tappetino ottimale per la rispettiva applicazione prima di usare l'agitatore.

Usare la Tabella 53 per determinare l'idoneità del materiale da laboratorio per l'utilizzo con i tappetini antiscivolo.
Prodotti di plastica / vetro	Indice per il potere aderente	Metallo	Indice per il potere aderente
Vetro (vetro borosilicato)	5	Acciaio inossidabile	5
Polipropilene (PP)	3	Alluminio	5
Polipropilene copolimero (PPCO)	3	Alluminio anodizzato	4
Polietilene (HPDE) (LPDE)	3		
Policarbonato (PC)	4		
Polietilene tereftalato (PET)	4		
Polietilene tereftalato copoliestere (PETG)	4		
Polimetilpentene (PMP)	3		
Politetrafluoroetilene (PTFE)	0*		
Copolimero di tetrafluoroetilene e esafluoropropilene (FEP)	0*		
Polimeri perfluoroalcossili (PFA)	2		
Silicone	1*		
Indici: 0 = nessun potere ad con indice 0 o 1 (PTFE, FEP	erente; 5 = massimo ? e silicone).	ootere aderente / *	Non usare recipienti

Tabella 53: Compatibilità dei materiali dei recipienti

AVVISO Tenere presente che la massima velocità di agitazione può essere raggiunta solo con materiali con indice «5» nella Tabella 53.

AVVISO Per materiali diversi dal vetro l'utente deve provare e determinare da sé quale tipo di materiale può essere usato per quale carico, velocità, durata di funzionamento e temperatura.

Campo di velocità e limite della durata di funzionamento per recipienti di vetro (riempiti al 30 %)

Valori limite per la velocità e la durata di funzionamento a temperatura ambiente per recipienti di vetro (riempiti al 30 %). Temperature superiori riducono il potere aderente. Un cambiamento della temperatura può comportare il distacco dei recipienti dal tappetino. Iniziare il processo di agitazione con recipienti acclimatati.

	Micropiastre / DeepWell	Becher			
	96 x 2 ml	25 ml	125 ml	250 ml	500 ml
125 giri/min	nessun limite di durata	1			
150 giri/min	72 ore	72 ore	72 ore	72 ore	72 ore
200 giri/min	72 ore	72 ore	72 ore	72 ore	72 ore
250 giri/min	72 ore				
300 giri/min					
350 giri/min					
400 giri/min					

 Tabella 54: Tappetino a bassa adesione - Campo di velocità e limite di durata a temperatura

 ambiente

	Micropiastre / DeepWell	Erlenm	neyer						
	96 x 2 ml	25 ml	125 ml	250 ml	500 ml	1000 ml	2000 ml	3000 ml	5000 ml
125 giri/ min	nessun limite di d	durata							
150 giri/ min	72 ore	72 ore	72 ore	72 ore	72 ore	72 ore	72 ore	72 ore	72 ore
200 giri/ min	72 ore	72 ore	72 ore	72 ore	72 ore	72 ore	72 ore	72 ore	72 ore
250 giri/ min	72 ore	72 ore	72 ore	72 ore	72 ore	72 ore	72 ore	48 ore	24 ore
300 giri/ min	72 ore	72 ore	72 ore	72 ore	72 ore	72 ore	48 ore	24 ore	
350 giri/ min	48 ore	48 ore	48 ore	48 ore	48 ore	48 ore	24 ore		
400 giri/ min	24 ore	24 ore	24 ore	24 ore	24 ore				

 Tabella 55: Tappetino ad alta adesione - Campo di velocità e limite di durata a temperatura

 ambiente

Preparazione

ATTENZIONE Per la manipolazione di prodotti di vetro indossare sempre occhiali e guanti protettivi.

Prima di posizionare i recipienti sul tappetino antiscivolo:

- 1. Controllare se i recipienti presentano danni. Non usare mai recipienti danneggiati.
- 2. Asciugare il recipiente con un panno pulito e asciutto per rimuovere tutta l'umidità.
- 3. Rimuovere ogni traccia di grasso e di altre impurità.
- Controllare che il fondo del recipiente e il tappetino antiscivolo siano asciutti e puliti.

ATTENZIONE Anche un film quasi non visibile o una singola goccia di acqua sul fondo del recipiente possono comportare che il recipiente perda l'adesione e si stacchi dal tappetino dopo un determinato intervallo di funzionamento.



Figura 97: Tappetino antiscivolo: asciugare il recipiente

5. Ridurre il volume di riempimento o chiudere i recipienti per evitare la fuoriuscita di liquidi.

ATTENZIONE Anche se il recipiente aderisce perfettamente, già una sola goccia di liquido che scorre giù sulla parete esterna può causare il distacco del recipiente dal tappetino dopo un certo periodo di funzionamento.

Posizionamento dei recipienti

- 1. Posizionare i recipienti sul tappetino e premerli verso il basso.
- 2. Spingere o tirare con delicatezza per assicurarsi che i singoli recipienti aderiscano affidabilmente al tappetino.

AVVISO Tenere presente che a temperature basse oppure durante un funzionamento prolungato può formarsi acqua di condensa in grado di causare il distacco accidentale dei recipienti. La temperatura del campione e del recipiente non devono differire dal campo di temperatura specificato per l'agitatore. Per informazioni dettagliate sul possibile campo di temperatura dell'agitatore utilizzato consultare «1. 1. Dati tecnici» a pagina 12.

Rimozione dei recipienti

- 1. Premere il pulsante Arresto sul display oppure attendere la fine del ciclo.
- 2. Attendere l'arresto completo della piattaforma.
- Tirare il collo del recipiente con delicatezza e uniformemente verso il lato e aspettare alcuni secondi. I recipienti con una superfici di fondo grande potrebbero metterci qualche secondo prima di staccarsi.



Figura 98: Tappetino antiscivolo: spingere o tirare con cautela per staccare il recipiente

ATTENZIONE Spingere o tirare con delicatezza senza esercitare grande forza fisica! I recipienti potrebbero staccarsi all'improvviso e versare il loro contenuto!

4. Se non è possibile staccare un recipiente, versare un po' di acqua intorno al fondo del recipiente.

Fare asciugare il tappetino all'aria prima di riutilizzarlo.

Rimozione del tappetino

1. Per staccare il tappetino antiscivolo sollevarlo su un lato e tirarlo lentamente, staccandolo dalla piattaforma, come illustrato nella Figura 99.



Figura 99: Tappetino antiscivolo: sollevare e tirare via il tappetino per rimuoverlo

Pulizia e disinfezione

Se il potere aderente diminuisce pulire il tappetino.

ATTENZIONE Non usare solventi, agenti detergenti o di lavaggio diversi da sapone delicato per piatti senza additivi per la cura della pelle. La concentrazione raccomandata è di massimo 2 ml di detersivo su 1 litro di acqua.

Accessori

AVVISO Il contatto diretto del tappetino con il detersivo concentrato riduce il potere aderente. In questo caso non è sufficiente sciacquare soltanto il tappetino. Se viene usato direttamente il detersivo, sciacquare il tappetino, metterlo per 24 ore ammollo in acqua e quindi farlo asciugare per 24 ore per ripristinare l'intero potere di adesione.

AVVISO Disinfettare solo con isopropanolo al 70 % o con etanolo al 70 %.

Leggermente sporco o impolverato:

- 1. Passare un panno umido o disinfettare il lato superiore del tappetino posato sulla piattaforma.
- 2. Inumidire il tappetino.
- 3. Eliminare i residui con una spugna.
- 4. Passare una spugna imbevuta di acqua pulita.
- 5. Disinfettare, se necessario.
- 6. Fare asciugare per 24 ore.

Fortemente sporco:

- 1. Rimuovere il tappetino dalla piattaforma per effettuare una pulizia completa.
- Mettere il tappetino ammollo in acqua addizionata di un detersivo diluito per piatti.
- 3. Sfregare accuratamente il tappetino con una spugna per rimuovere i residui.
- 4. Dopo la pulizia, sciacquare il tappetino abbondantemente con acqua.
- 5. Disinfettare, se necessario.
- 6. Fare asciugare per 24 ore.
- Posizionare il tappetino come spiegato alla sezione «Posizionamento del tappetino asciutto» a pagina 108 oppure «Posizionamento del tappetino bagnato» a pagina 108.

3. 3. 12. Installazione del collettore del gas

Il collettore del gas permette l'iniezione di gas inerte, come ad es. azoto o anidride carbonica, in diversi recipienti durante il funzionamento dell'agitatore (vedere il capitolo «1. 2. 12. Accessori generali» a pagina 43 per le informazioni di ordinazione).

Il collettore del gas è un'opzione installata in fabbrica per gli agitatori delle serie Solaris 2000 I/R e 4000 I/R. Questo collettore è anche disponibile come accessorio da installare successivamente dall'utente se non installato originariamente sul dispositivo. Tutti gli agitatori Solaris 2000 I/R e 4000 I/R hanno un foro passacavo su una delle due pareti laterali (vedere il capitolo «1. 4. Informazioni generali sul prodotto» a pagina 45 per informazioni sulla posizione) destinato al passaggio del tubo di alimentazione del gas da laboratorio al gruppo collettore. Il collettore del gas possiede un raccordo di ingresso per l'alimentazione del gas e otto raccordi di uscita per i recipienti.



- Tubo di alimentazione del gas da laboratorio
- ② Gruppo collettore del gas
- ③ Raccordi di uscita con tubi flessibili
- ④ Raccordo di ingresso
- 5 Viti di fissaggio

Figura 100: Collettore del gas installato sull'agitatore Solaris 4000 R

Per l'installazione procedere come segue:

- 1. Fissare accuratamente il collettore del gas ② sulla parete posteriore della camera dell'agitatore con le viti ⑤ fornite.
- Rimuovere il tappo di plastica dal lato esterno della boccola del passacavo (vedere il capitolo «1. 4. Informazioni generali sul prodotto» a pagina 45 per informazioni sulla posizione).
- 3. Con una taglierina, praticare un'incisione a croce nella membrana della boccola di gomma all'interno della camera dell'agitatore.

- 4. Tagliare un pezzo di tubo flessibile alla lunghezza adatta per usarlo come tubo di alimentazione del gas da laboratorio ①.
- 5. Collegare il tubo di alimentazione del gas da laboratorio ① al raccordo di ingresso scanalato ④ sul collettore.
- Assicurarsi che il tubo di alimentazione del gas da laboratorio ① stia a distanza dalle leve della cappa e farlo passare attraverso la boccola di gomma nel foro passacavo verso l'esterno.
- 7. Collegare il tubo di alimentazione del gas da laboratorio ① alla fonte del gas.
- Tagliare 8 pezzi di tubo flessibile e collegarli alle uscite 3 del collettore. Per un numero di recipienti superiore a 8 utilizzare connettori a «Y»; per un numero di recipienti inferiore a 8, chiudere a morsetto i tubi non utilizzati per ridurre la perdita di gas.
- 9. Usare tappi e guarnizioni adatti per inserire i tubi nei recipienti.

3.4. Caricamento e uso previsto



Rischio di incendio causato da reazioni chimiche innescate. Non usare l'agitatore a velocità che provocano la fuoriuscita dei contenuti dai vassoi.

AVVERTENZA

Aumentare lentamente la velocità. Provare con l'acqua prima di usare le sostanze chimiche.



La sicurezza può essere pregiudicata da un caricamento non corretto e da accessori difettosi.

ATTENZIONE

- Assicurarsi che il carico (accessori e campioni) siano collocati in maniera simmetrica rispetto al centro della piattaforma. Con l'uso di una piattaforma a due livelli, assicurarsi che la piattaforma inferiore abbia un carico maggiore rispetto alla piattaforma superiore.
- Non sovraccaricare mai l'agitatore. Per il carico massimo consultare «Dati tecnici» a pagina 12. Il carico comprende il peso della piattaforma montata sull'agitatore, degli accessori e dei campioni.
- Assicurarsi che gli accessori siano montati correttamente prima di usare l'agitatore. Attenersi alle istruzioni riportate al capitolo «Accessori» a pagina 89.



Rischio di lesioni alle dita o alla mano causate dalla cappa che cade giù.

- ATTENZIONE
- La cappa dell'agitatore potrebbe cadere giù e chiudersi all'improvviso e schiacciare o incastrare le dita o le mani, provocando lesioni. Ciò può accadere se la molla della cappa è difettosa e non riesce a mantenere la cappa nella posizione aperta.

Non utilizzare l'agitatore con una molla della cappa difettosa. Chiamare il servizio di assistenza della Thermo Fisher Scientific e fare sostituire subito la molla della cappa.

Caricamento

Assicurarsi che il carico, compresi gli accessori e i campioni, sia collocato in maniera simmetrica rispetto al centro della piattaforma. Quando viene usata una piattaforma a due livelli senza sfruttare l'intera capacità, tenere presente quanto segue: per ottenere la migliore performance possibile, caricare preferibilmente la piastra inferiore e disporre il carico in maniera simmetrica rispetto al centro della piattaforma.

Con l'uso di un agitatore a termoregolazione assicurarsi di posizionare il carico in modo che alla chiusura della cappa e con la cappa chiusa vi sia uno spazio libero sufficiente per evitare collisioni e possibili danni al carico.



Figura 101: Esempio di una piattaforma caricata bene

												ā.	- A
•					*			٠	٠	÷	-	•	
•	٠	٠	٠	٠	٠	٠	٠		٠		. "	÷	·(•)•
		٠	*	٠	*	*	*	*	*	٠	٠	*	
			٠	*		*		*	*		×		· 🙈 📢 🕩
٠		٠	*	•	*	*	*	٠	*	°,	٠	*	\bigcirc
	×						٠	•			*		
		•			*		*	٠	٠				· 👝 🜔 🕽
•						*	٠			*	٠		
٠								*			٠	٠	
•		•			*		*		*	*	٠	*	$\cdot \land (\cdot)$
•	٠	•	•	•	•	•	•	٠	•	٠		٠	

Figura 102: Esempio di una piattaforma caricata male

AVVISO II massimo carico comprende il peso della piattaforma montata sull'agitatore, degli accessori e dei campioni. Per maggiori informazioni sul massimo carico vedere «1. 1. Dati tecnici» a pagina 12.

Uso previsto

Le informazioni riguardanti i limiti di velocità-carico sono solo valori indicativi per l'utilizzo di questo prodotto. Un funzionamento uniforme con un basso livello di vibrazioni o in assenza di vibrazioni dipende da molteplici fattori, quali ad esempio il livello di riempimento, il tipo e le condizioni dei supporti e dei tappetini, il tipo di recipiente, la sistemazione dei recipienti sulla piattaforma, la superficie di appoggio dell'agitatore e la velocità impostata.

Quando viene impostata una combinazione nuova o sconosciuta oppure se viene modificato uno di questi parametri, tenere presente quanto segue: aumentare lentamente la velocità sopra 175 giri/min. e controllare se si presentano vibrazioni indesiderate del carico o movimenti indesiderati dell'agitatore. Il cliente stesso è responsabile di un funzionamento sicuro dell'agitatore.

Gli agitatori della serie Solaris sono progettati appositamente in modo da emettere un basso livello di calore e di vibrazioni, il che li rende particolarmente idonei per l'impiego in una serie di dispositivi da laboratorio, compresi camere climatiche, incubatori e refrigeratori da laboratorio. In considerazione delle svariate condizioni ambientali specifiche, come ad esempio la stabilità dei ripiani e supporti, in caso dell'impiego all'interno di dispositivi da laboratorio l'utente è responsabile di un funzionamento sicuro. Aumentare lentamente la velocità e fare attenzione a movimenti indesiderati del dispositivo.

Velocità sicura

Per gli agitatori della serie Solaris la velocità sicura indipendente dal carico è riportata nella Tabella 56. Il carico comprende il peso della piattaforma montata sull'agitatore, degli accessori e dei campioni. Per maggiori informazioni sul carico massimo consultare «Dati tecnici» a pagina 12.

Тіро	Velocità max. sicura
Solaris 2000	175 giri/min
Solaris 4000	175 giri/min
Solaris 2000 I	250 giri/min
Solaris 2000 R	250 giri/min
Solaris 4000 I	250 giri/min
Solaris 4000 R	250 giri/min

Tabella 56: Velocità massima sicura

Velocità massima

Alla velocità massima di 525 giri/min. gli agitatori della serie Solaris possono essere usati con un carico limitato secondo la Tabella 57. Il carico comprende il peso della piattaforma montata sull'agitatore, degli accessori e dei campioni. Tenere conto delle velocità ammesse per i supporti e recipienti. Questi potrebbero non essere compatibili con questa velocità.

Тіро	Carichi max. sicuri
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

Tabella 57: Carico massimo alla velocità massima

AVVISO Tenere sempre presente che con l'aumento della velocità i supporti potranno iniziare ad aprirsi provocando temporaneamente vibrazioni supplementari del sistema a seguito di un maggiore movimento dei recipienti.

Grafici peso/velocità

I grafici peso/velocità riportati di seguito aiutano a impostare la combinazione velocità-carico ottimale per l'uso previsto. L'area verde mostra le combinazioni di velocità-carico che non comportano vibrazioni o causano vibrazioni minime durante l'agitazione. L'area rossa mostra le combinazioni di velocità-carico che potrebbero causare forti vibrazioni durante l'agitazione e che potrebbero provocare uno spostamento indesiderato dell'agitatore. Questi diagrammi sono forniti a titolo puramente indicativo e pertanto l'operatore deve fare attenzione quando la combinazione di velocità-carico si avvicina all'area rossa.

Tenere presente che le condizioni dell'applicazione specifica potrebbero causare un comportamento indesiderato del carico o del dispositivo prima di raggiungere la linea di confine mostrata. Aumentare gradualmente la velocità per valutare il comportamento del carico specifico. Il carico comprende il peso della piattaforma montata sull'agitatore, degli accessori e dei campioni.

AVVISO Ogni linea blu orizzontale nei grafici in basso indica il peso a vuoto di una piattaforma specifica.

Solaris 2000



Figura 103: Solaris 2000 – Uso previsto



Solaris 4000

Figura 104: Solaris 4000 – Uso previsto

Caricamento e uso previsto



Figura 105: Solaris 2000 I / 2000 R - Uso previsto



Solaris 4000 I / 4000 R

Figura 106: Solaris 4000 I / 4000 R - Uso previsto

4. Manutenzione e cura



AVVERTENZA

Rischi provocati dalla manipolazione di sostanze pericolose

- Per l'agitazione di materiali pericolosi attenersi al manuale di biosicurezza nei laboratori («Laboratory Biosafety Manual») dell'Organizzazione Mondiale della Sanità (OMS) e alle norme nazionali vigenti in materia. Se vengono agitate prove microbiologiche del gruppo di rischio II (secondo il manuale «Laboratory Biosafety Manual» dell'organizzazione di salute mondiale WHO) devono essere impiegate chiusure bio-aerosol con tenuta aerosol. Sul sito dell'Organizzazione Mondiale della Sanità OMS (www.who.int); (WHO, World Health Organisation) cercare il manuale «Laboratory Biosafety Manual». Con materiali con un gruppo di rischio più alto deve essere previsto più di un provvedimento di protezione.
- Se l'agitatore o i suoi componenti sono stati contaminati da sostanze tossiche o patogene, è necessario adottare misure di disinfezione idonee («Decontaminazione» a pagina 125; «Disinfezione» a pagina 125).
- Al verificarsi di una situazione di pericolo, disinserire l'alimentazione dell'agitatore e abbandonare subito l'ambiente circostante l'agitatore.



Danni alla salute provocati da sostanze infettive

AVVERTENZA

Qualora liquidi o altri materiali dovessero essere versati accidentalmente sotto la piattaforma, spegnere subito l'agitatore, staccare la spina e rimuovere la piattaforma («Piattaforme» a pagina 19).

Eliminare il materiale versato seguendo le regolari procedure adottate nel laboratorio. Utilizzare i dispositivi di protezione individuale idonei.

Tutte le regolazioni interne e le riparazioni devono essere effettuate da un tecnico autorizzato del servizio di assistenza. Il corpo dell'agitatore non deve essere aperto dall'operatore.

Attenersi a tutta la documentazione fornita insieme al rispettivo accessorio e riportante le specifiche informazioni sulla manutenzione e sulla pulizia accurate. Usare le informazioni contenute nel presente capitolo solo a titolo indicativo.

4.1. Principi fondamentali

Per la protezione di persone, ambiente e materiali è obbligatorio pulire e, se necessario, disinfettare periodicamente l'agitatore e gli accessori.

Thermo Fisher Scientific raccomanda di pulire e disinfettare manualmente l'agitatore da laboratorio almeno una volta al mese. La normale aria nell'ambiente chiuso contiene migliaia di microrganismi circolanti in grado di insediarsi nell'agitatore e di mettere a rischio le colture.

Thermo Fisher Scientific raccomanda l'uso di etanolo al 70 % o di isopropanolo al 70 % o di un disinfettante a base di sali quaternari di ammonio al 10 % o meno.

AVVISO Il meccanismo può essere danneggiato da liquidi penetranti. Badare affinché i liquidi, in particolare i solventi organici, non vengano a contatto con il meccanismo o con il cuscinetto del meccanismo. I solventi organici alterano il grasso dei cuscinetti del meccanismo. In caso di liquidi che penetrano in componenti dell'agitatore non accessibili per la pulizia, non usare l'agitatore e contattare il servizio di assistenza post vendita.

AVVISO Procedure o prodotti non omologati possono aggredire i materiali dell'agitatore e provocare malfunzionamenti. Non utilizzare procedure di pulizia o decontaminazione diverse, se non si è sicuri che queste siano idonee per l'apparecchiatura. Utilizzare solo detergenti che non danneggiano l'apparecchiatura. In caso di dubbio rivolgersi al produttore del detergente.

AVVISO Sugli agitatori a termoregolazione assicurarsi di lasciare aperta la cappa per fare evaporare gli agenti detergenti e l'acqua.

- Staccare la spina di alimentazione elettrica prima di effettuare la pulizia, disinfezione o decontaminazione.
- Rimuovere dall'agitatore gli accessori e la piattaforma montati prima di effettuare la pulizia, disinfezione o decontaminazione.
- Usare acqua calda con un detergente delicato e un panno morbido per pulire i materiali. In caso di dubbio rivolgersi a Thermo Fisher Scientific. Sciacquare con acqua pulita ed asciugare accuratamente.
- Non utilizzare mai detergenti corrosivi, come acido fosforico, candeggina o polvere abrasiva.
- Utilizzare solo disinfettanti con un valore pH di 6–8.
- Eliminare subito eventuali liquidi traboccati con un panno che non sfilaccia inumidito con un detergente non corrosivo (vedere le specifiche del produttore del detergente).
- Liquidi versati possono penetrare sotto la piattaforma. Se un liquido traboccato scorre sotto la piattaforma, smontare la piattaforma ed eliminare accuratamente il liquido dalla vaschetta sotto la piattaforma. Disinfettare o decontaminare, se necessario.
- Controllare i componenti dell'agitatore ed eliminare eventuali impurità e liquidi traboccati.

Ispezione degli accessori

AVVISO Non utilizzare un agitatore o accessori che presentano segni di danneggiamento. Si raccomanda di ispezionare gli accessori periodicamente nell'ambito della routine di manutenzione per garantire la sicurezza.

Dopo un'accurata pulizia degli accessori questi devono essere controllati per rilevare eventuali danni, segni di usura e corrosione.

Componenti di metallo

In caso di danni, come ad esempio corrosione, usura o incrinature, gli accessori devono essere messi fuori servizio subito.

Componenti in plastica

Controllare se vi sono segni di incrinature, usura, graffi e crepe. In caso di danni, l'oggetto ispezionato deve essere messo fuori servizio subito.

Tappetini antiscivolo

Prima di ogni utilizzo ispezionare i tappetini antiscivolo per individuare eventuali segni di usura o danni, compresi screpolature, scoloriture, ingiallimenti, infragilimento, deformazioni, abrasioni superficiali o attacchi chimici. Scartare subito i tappetini antiscivolo che presentano segni di usura o di danneggiamento.

4.2. Pulizia

AVVISO Prima di adottare una procedura di pulizia, accertarsi presso il produttore del detergente che la procedura prevista non danneggi l'apparecchiatura.

AVVISO II meccanismo può essere danneggiato da liquidi penetranti. Badare affinché i liquidi, in particolare i solventi organici, non vadano a finire sul meccanismo o sul cuscinetto del meccanismo. I solventi organici alterano il grasso dei cuscinetti del meccanismo. Lavare la parte esterna del dispositivo con un panno morbido e una soluzione di acqua e sapone neutro, sciacquare con acqua pulita e asciugare accuratamente.

Per informazioni sulla pulizia accurata dell'agitatore e degli accessori utilizzati consultare «Principi fondamentali» a pagina 123.

AVVISO Sugli agitatori a termoregolazione assicurarsi di lasciare aperta la cappa per fare evaporare gli agenti detergenti e l'acqua.

Touch-screen

Per la pulizia del touch-screen:

- 1. Staccare la spina di alimentazione elettrica.
- 2. Pulire lo schermo touch-screen con un panno in microfibra asciutto.
- 3. Se necessario, inumidire il panno in microfibra con acqua e passarlo un'altra volta sul touch-screen.
- 4. Lasciare asciugare all'aria.

4.3. Disinfezione

AVVERTENZA Rischi provocati dalla manipolazione di sostanze pericolose. Non toccare parti contaminate. Pericolo di infezione al contatto con i componenti contaminati. In caso di rottura del recipiente o versamenti del contenuto, materiale infettivo può penetrare nell'agitatore. In caso di contaminazione assicurarsi che non venga messo in pericolo nessuno. Decontaminare subito tutte le parti coinvolte.

AVVISO Danneggiamento delle apparecchiature causato da detergenti o metodi di disinfezione non idonei. Assicurarsi che il metodo o l'agente disinfettante non danneggino l'apparecchiatura. In caso di dubbio rivolgersi al produttore del disinfettante. Attenersi alle avvertenze di sicurezza e alle istruzioni per la manipolazione dei disinfettanti utilizzati.

- 1. Passare su tutti i componenti e su tutte le superfici etanolo al 70 % come richiesto dal rispettivo livello di disinfezione. Non bagnare le aree con componenti elettronici esposti.
- 2. Lasciare asciugare all'aria.

AVVISO Sugli agitatori a termoregolazione assicurarsi di lasciare aperta la cappa per fare evaporare gli agenti detergenti e l'acqua.

L'operatore stesso è responsabile che venga raggiunto il livello di disinfezione richiesto dagli standard specifici.

4.4. Decontaminazione

AVVERTENZA Rischi provocati dalla manipolazione di sostanze pericolose. Non toccare parti contaminate. Pericolo di contaminazione al contatto con i componenti contaminati. In caso di rottura della provetta o versamenti del contenuto, materiale contaminato può penetrare nell'agitatore. In caso di contaminazione assicurarsi che non venga messo in pericolo nessuno. Decontaminare subito tutte le parti coinvolte.

AVVISO Danneggiamento delle apparecchiature causato da detergenti o metodi di decontaminazione non idonei. Assicurarsi che il metodo o l'agente di decontaminazione non danneggino l'apparecchiatura. In caso di dubbio rivolgersi al produttore dell'agente di decontaminazione. Attenersi alle avvertenze di sicurezza e alle istruzioni per la manipolazione degli agenti decontaminanti utilizzati.

Il seguente metodo è consigliato da Thermo Fisher Scientific.

- 1. Passare etanolo al 70 % su tutti i componenti e tutte le superfici. Non bagnare le aree con componenti elettronici esposti.
- 2. Lasciare asciugare all'aria.

AVVISO Sugli agitatori a termoregolazione assicurarsi di lasciare aperta la cappa per fare evaporare gli agenti detergenti e l'acqua.

L'operatore stesso è responsabile che venga raggiunto il livello di decontaminazione richiesto dagli standard specifici.

4.5. Autoclavaggio

AVVISO Non superare mai i valori ammessi per quanto riguarda la temperatura e la durata dell'autoclavaggio. Non sono ammessi additivi chimici nel vapore.

Prima dell'autoclavaggio:

- Disassemblare tutti gli accessori.
- Sciacquare accuratamente per eliminare ogni traccia di sostanze chimiche o detergenti. Effettuare un risciacquo finale con acqua distillata.

Per informazioni riguardanti l'autoclavabilità consultare la seguente tabella:

Pos.	Autoclava- bile	Specifiche
Agitatore	No	-
Piattaforma	Sì	121 °C, 20 min
Supporto porta beute (senza bussole a molla)	Sì	121 °C, 20 min
Cacciavite	No	-
Supporto per micropiastra / piastra DeepWell	Sì	121 °C, 20 min
Tappeto antiscivolo	No	-
Rack porta becher	Sì	121 °C, 20 min
Supporto per imbuto separatore	Sì	121 °C, 20 min
Rack portaprovette - componenti di metallo	Sì	121 °C, 20 min
Rack portaprovette - componenti di plastica	No	-
Rack portaprovette – inserti in materiale espanso	No	-
Supporto regolabile per recipienti - parti di metallo	Sì	121 °C, 20 min
Supporto regolabile per recipienti - tappetino antiscivolo	No	-
Supporto multiuso - supporto in metallo	Sì	121 °C, 20 min
Supporto multiuso - tappetino di gomma	No	-
Supporto ad angolo regolabile per rack portaprovette -	Sì	121 °C, 20 min
componenti di metallo		
Supporto ad angolo regolabile per rack portaprovette -	No	-
componenti di plastica		
Supporto ad angolo regolabile per rack portaprovette -	No	-
inserti in materiale espanso		

Tabella 58: Autoclavabilità dei materiali

Assicurarsi che venga raggiunta la sterilità necessaria in conformità ai requisiti specifici.

L'aspetto e il colore possono cambiare leggermente dopo l'autoclavaggio.

4.6. Calibrazione di temperatura

Il valore di temperatura nominale di un agitatore a termoregolazione può essere calibrato con un dispositivo di misurazione della temperatura di precisione nota per garantire la piena riproducibilità del processo di agitazione (incubato) o refrigerato (raffreddato e riscaldato).

4. 6. 1. Buone pratiche per la calibrazione della temperatura

Seguire accuratamente le istruzioni prima e durante la calibrazione della temperatura sull'agitatore. Qualsiasi disturbo o deviazione dalle buone pratiche raccomandate può impedire lo stabilizzarsi della temperatura del campione e causare il fallire del ciclo di calibrazione.

Le buone pratiche prevedono di rispettare quanto segue:

- Usare un matraccio con un volume di 250 ml.
- Riempire il matraccio esattamente con 100 ml di acqua.
- Chiudere la cappa completamente e mantenerla chiusa per l'intero ciclo di calibrazione.
- Fare passare un intervallo di tempo sufficiente per permettere che la temperatura del campione si sia stabilizzata completamente.
- Fare attenzione che le condizioni ambientali, come la temperatura o l'umidità, non cambino durante il ciclo di calibrazione. Perfino cambiamenti piccoli, come l'apertura di una porta o di una finestra o variazioni alle impostazioni del sistema di condizionamento aria, possono pregiudicare il risultato del processo di calibrazione in corso.
- Si raccomanda di calibrare l'agitatore una volta all'anno.

4. 6. 2. Effettuazione della calibrazione di temperatura

Procedere come segue per effettuare la calibrazione:

- 1. La calibrazione può essere avviata in tre modi diversi:
 - » Alla fine del processo di configurazione iniziale descritto in «Configurazione iniziale» a pagina 55 si ha la possibilità di avviare la calibrazione della temperatura invece di uscire dal processo di configurazione iniziale.
 - » L'agitatore visualizza un promemoria per la calibrazione annuale.
 - » Quindi la calibrazione può essere avviata toccando il simbolo Impostazioni nella barra dei menu, seguito dal pulsante Calibrazione più in fondo sulla schermata Calibrazione temp che appare.
 - » Si può toccare il simbolo **Impostazioni** nella barra dei menu, seguito dal pulsante **Calibrazione** per avviare il processo in qualsiasi momento ciò risulti necessario.

AVVISO Il promemoria di calibrazione può essere disattivato spuntando la casella **Disattiva avvisi di calibrazione** sulla schermata **Impostazioni**.

2. Attendere fino a quando appare la schermata **Calibrazione temp** illustrata nella Figura 107.



Figura 107: Calibrazione: schermata iniziale

3. Toccare il pulsante Avvia calibrazione per iniziare la procedura.



Figura 108: Calibrazione: richiesta di riempire un matraccio di Erlenmeyer

- 4. Seguire le istruzioni visualizzate sulla schermata: Riempire il matraccio di Erlenmeyer da 250 ml con 100 ml di acqua.
- 5. Posizionare il matraccio in un supporto e fissare il supporto all'incirca al centro della piattaforma di agitazione.
- Appendere la sonda di un dispositivo di misurazione della temperatura con accuratezza nota nel matraccio. La sonda deve essere immersa nel liquido, senza toccare il fondo o i lati del matraccio.
- 7. Chiudere il matraccio per evitare la fuoriuscita del liquido.
- 8. Chiudere la cappa sopra la piattaforma di agitazione.

9. Toccare il pulsante Avanti per continuare.



Figura 109: Calibrazione: richiesta di avviare la calibrazione

AVVISO Appena viene toccato il pulsante Avvio l'agitatore inizia subito con l'agitazione.

 Seguire le istruzioni visualizzate sulla schermata: Toccare il pulsante Avvio per usare la temperatura preimpostata oppure toccare il campo Valore nominale temperatura e inserire la temperatura desiderata prima di toccare il pulsante Avvio.

La schermata visualizza come la camera dell'agitatore viene riscaldata (o raffreddata) alla temperatura di calibrazione selezionata.



Figura 110: Calibrazione: Stabilizzazione della camera alla temperatura di calibrazione

11. Attendere almeno due ore e mezza (150 minuti) finché la misurazione della temperatura nella camera non abbia raggiunto condizioni stabili.

Le condizioni stabili sono indicate dal messaggio La temperatura nella camera si è stabilizzata, come illustrato nella Figura 111 in basso.



Figura 111: Calibrazione: Camera stabilizzata con successo alla temperatura di calibrazione

- 12. Seguire le istruzioni visualizzate sulla schermata:
 - » Appena viene visualizzata la schermata illustrata nella Figura 111, confrontare a intervalli regolari la temperatura rilevata dalla termocoppia. Se ritenete che la temperatura sia stabile si può o toccare il pulsante Modifica per correggere la temperatura oppure toccare il pulsante Fine per terminare il processo di calibrazione.



Figura 112: Calibrazione: Correzione della temperatura visualizzata dall'agitatore

- Per mezzo della rotella di selezione che appare correggere la temperatura visualizzata dall'agitatore, inserendo il valore rilevato dal dispositivo di misurazione di precisione.
- 14. Toccare il pulsante **Salva** per salvare le modifiche apportate.

15. Il messaggio **Calibrazione temperatura completata** indica che il processo è stato completato con successo.



Figura 113: Calibrazione: Calibrazione della temperatura completata con successo

Si può toccare Ricalibra per ripetere la procedura di calibrazione.

4. 6. 3. Individuazione della temperatura calibrata

Si ha la possibilità di consultare in qualsiasi momento l'attuale temperatura calibrata:

- 1. Toccare il simbolo **Impostazioni** nella barra dei menu, seguito dal pulsante **Calibrazione** più in fondo sulla schermata **Calibrazione temp** che appare.
- 2. Appare la schermata **Calibrazione temp** illustrata nella Figura 107. Questa schermata visualizza l'attuale **Temperatura agitatore calibrata**.

AVVISO In alternativa, gli agitatori a termoregolazione permettono di consultare la temperatura di calibrazione sulla schermata di impostazione della temperatura. Un esempio viene illustrato nella sezione «Impostazione della temperatura» a pagina 65.

4.7. Installazione del firmware

Toccando il pulsante Firmware sulla schermata File & Info (vedere «File & Info» a pagina 78) il sistema chiede all'utente di inserire il codice da amministratore. Dopo l'inserimento del codice di accesso si apre la schermata Firmware illustrata nella Figura 114. Questa schermata permette di installare un firmware nuovo da una chiavetta USB.

In base agli accordi di assistenza stipulati è possibile che si riceva dal tecnico di assistenza post vendita un link per scaricare il nuovo firmware. Fare attenzione a installare solo un firmware proveniente da una fonte autorizzata. Un software proveniente da una fonte non autorizzata può comportare l'estinzione della garanzia concessa da Thermo Fisher Scientific.

- 1. Scaricare il file zip dal link fornito dal proprio tecnico di assistenza post vendita.
- 2. Decomprimere il file zip su una chiavetta USB vuota.

AVVISO Accertarsi che la chiavetta USB disponga di uno spazio di memoria sufficiente per consentire l'estrazione del file zip.

- Se l'agitatore è ancora in funzione, arrestarlo, premendo il pulsante Arresto sul display.
- 4. Confermare eventuali allarmi presenti, come descritto alla sezione «Stato» a pagina 67.

AVVISO Se si cerca di effettuare un aggiornamento mentre l'agitatore è in funzione, il sistema emette un messaggio per informare l'utente che Non è permesso eseguire un aggiornamento del firmware mentre l'agitatore si trova in una fase di agitazione attiva. Questo messaggio ha un pulsante di Arresta Shaker per interrompere il ciclo ed eseguire l'aggiornamento.

- 5. Toccare il simbolo File & Info nella barra di navigazione.
- 6. Scorrere fino in fondo alla schermata **File & Info** per vedere il pulsante **Firmware**.
- 7. Toccare il pulsante **Firmware** per accedere alla schermata Firmware che visualizza le versioni firmware attualmente installate.
- Inserire il dispositivo USB in una delle porte USB dell'agitatore. Per individuare la posizione delle porte USB consultare la sezione «Informazioni generali sul prodotto» a pagina 45.
- 9. Inserire il codice di accesso da amministratore.



Figura 114: Schermata del firmware

10. Toccare il pulsante Aggiorna firmware.

AVVISO Se l'agitatore non trova nessuna chiavetta USB, il sistema chiaderà all'utente di Inserire una chiavetta USB con il firmware da installare per iniziare l'aggiornamento. Adesso inserire la chiavetta USB o controllare se è inserita correttamente nella porta USB e porre rimedio al problema. 11. Se è disponibile un firmware da aggiornare, sulla schermata Aggiornamento firmware appare la lista illustrata nella Figura 115. Toccare l'elemento firmware che si desidera aggiornare.

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

Figura 115: Selezione dell'elemento firmware da installare

AVVISO Gli elementi firmware che sono già aggiornati appaiono sbiaditi e non possono essere selezionati. Se nessuno degli elementi è da aggiornare, il sistema emette un messaggio per informare l'utente che Attualmente non sono presenti nuovi aggiornamenti per il software.

- 12. Toccare il pulsante Seleziona.
- Se la chiavetta di memoria contiene più di una versione firmware, il sistema chiede all'utente di selezionarne una. Toccare la versione che si desidera selezionare e quindi toccare Seleziona.





AVVISO In caso di dubbi sulla versione da installare, rivolgersi al tecnico di assistenza post vendita che ha fornito il link per il download.

14. Appare una lista degli elementi da aggiornare. Toccare **OK** per avviare l'aggiornamento.



Figura 117: Lista degli elementi da aggiornare

AVVISO L'interfaccia utente gestisce le dipendenze reciproche per l'utente: l'utente sceglie un elemento firmware per l'aggiornamento e il software aggiunge automaticamente tutti gli elementi necessari da installare.

AVVISO Se viene selezionata l'opzione H.M.I. dalla lista illustrata nella Figura 115 per aggiornare solo il firmware dell'interfaccia utente, per un periodo che può durare fino a un minuto dopo l'avvio dell'aggiornamento firmware il display dell'interfaccia utente non risponderà ai contatti sull'area attiva del touch-screen. <u>Non</u> spegnere l'agitatore durante questa fase. Altrimenti l'aggiornamento non può essere completato con successo. Dopo un breve intervallo di tempo l'agitatore viene riavviato e quindi è nuovamente pronto al funzionamento.

AVVISO Se vengono selezionate due o tre opzioni dalla lista illustrata nella Figura 115 per aggiornare <u>diversi</u> componenti firmware in una volta, l'agitatore <u>non deve</u> essere spento in nessun momento durante l'intero processo di aggiornamento.

15. Appaiono diversi messaggi che informano l'utente in merito al progresso dell'installazione, come illustrato nella Figura 118.



Figura 118: Messaggi durante l'installazione dell'aggiornamento firmware

16. Appena il sistema chiede all'utente di riavviare l'agitatore, spegnere e riaccendere l'agitatore per mezzo dell'interruttore generale.

4.8. Sostituzione della piattaforma

La procedura base per l'installazione di una piattaforma è descritta al capitolo «3. 3. 1. Installazione della piattaforma» a pagina 90. Una lista completa delle piattaforme di ricambio per i singoli agitatori si trova alla sezione «1. 2. 1. Piattaforme» a pagina 19.

AVVISO I modelli di agitatori a termoregolazione devono essere utilizzati solo con le piattaforme preinstallate in fabbrica.

4.9. Manutenzione

La Thermo Fisher Scientific consiglia di fare effettuare una volta all'anno una manutenzione dell'agitatore e degli accessori da parte del servizio di assistenza autorizzato. Il tecnico del servizio di assistenza controlla quanto segue:

- equipaggiamento elettrico
- l'idoneità del luogo di installazione
- sistema di sicurezza
- accessori utilizzati
- fissaggio dei supporti, delle piattaforme e degli accessori sull'agitatore

Prima di effettuare lavori di manutenzione e assistenza, l'agitatore e gli accessori devono essere puliti e decontaminati accuratamente per garantire un'ispezione completa e sicura.

Per queste prestazioni la Thermo Fisher Scientific offre contratti di ispezione e di assistenza tecnica. Le riparazioni eventualmente necessarie vengono effettuate gratuitamente nell'ambito del periodo di garanzia ed addebitate se fuori dalla garanzia. Questo vale solo, se gli interventi di manutenzione sull'agitatore sono stati eseguiti da un tecnico autorizzato del servizio di assistenza di Thermo Fisher Scientific.

4.10. Spedizione e smaltimento

AVVERTENZA Danni alla salute provocati da sostanze infettive. Se l'agitatore e gli accessori vengono messi fuori servizio per essere smaltiti, questi devono essere puliti e all'occorrenza disinfettati o decontaminati. In caso di dubbio contattare il servizio di assistenza della Thermo Fisher Scientific.

Per lo smaltimento dell'agitatore attenersi alle norme vigenti nel rispettivo Paese. Anche per domande riguardanti lo smaltimento dell'agitatore, contattare il servizio di assistenza Thermo Fisher Scientific. Le informazioni di contatto sono riportate sul retro delle presenti istruzioni oppure sul sito <u>www.thermofisher.com</u>.

Attenersi alle informazioni riguardanti il trasporto e la spedizione («Trasporto» a pagina 53, «Spedizione» a pagina 60).

5. Eliminazione guasti

N.	Descrizione	Soluzioni	Simbolo
Errori			
11999	Si è verificato un errore interno.	Si è verificato un errore interno. Riavviare il dispositivo, spegnendo e riaccendendo l'interruttore generale. Se il messaggio d'errore continua essere visualizzato, contattare un tecnico del servizio di assistenza.	
Avvertime	enti		
3	È richiesta la calibrazione della temperatura. È passato 1 anno.	Il periodo di validità di un anno è scaduto. Ripetere la calibrazione della temperatura come spiegato nella sezione «4. 6. Calibrazione di temperatura» a pagina 127.	CAL
4	Alimentazione elettrica ripristinata e avvio automatico abilitato.	Durante l'ultimo ciclo si è verificata un'interruzione dell'alimentazione. Dopo il ripristino dell'alimentazione elettrica il ciclo è stato continuato automaticamente.	Ŧ
6	Guasto ai sensori della temperatura ambiente.	La temperatura misurata dal sensore della temperatura ambiente non è plausibile.	III
10	Velocità per la ventola 1 supera il valore limite.	Velocità per la ventola 1 supera la soglia di plausibilità.	•
11	Velocità per la ventola 2 supera il valore limite.	Velocità per la ventola 2 supera la soglia di plausibilità.	•2
12	Velocità per la ventola 3 supera il valore limite.	Velocità per la ventola 3 supera la soglia di plausibilità.	•
13	Velocità per la ventola 4 supera il valore limite.	Velocità per la ventola 4 supera la soglia di plausibilità.	•2
20	Cappa è aperta da un periodo di tempo prolungato.	Chiudere la cappa dell'agitatore per mantenere la temperatura della camera.	
Allarmi			
1	Allarme di temperatura alta	La temperatura della camera ha superato il livello d'allarme superiore. Controllare i campioni, le condizioni ambientali e/o modificare le impostazioni.	III
2	Allarme di temperatura bassa	La temperatura della camera ha superato il livello d'allarme inferiore. Controllare i campioni, le condizioni ambientali e/o modificare le impostazioni.	
3	Il riavvio automatico non è riuscito dopo un'interruzione dell'alimentazione.	Durante l'ultimo ciclo si è verificata un'interruzione dell'alimentazione. Il riavvio automatico non è riuscito.	Ċ
5	Accelerazione azionamento troppo bassa. Non è possibile raggiungere la velocità nominale desiderata.	Non è stato possibile raggiungere la velocità impostata nel tempo previsto. Controllare le impostazioni e/o il carico (accessori e campioni) sulla piattaforma.	\odot

Spedizione e smaltimento

N.	Descrizione	Soluzioni	Simbolo
7	La misurazione della velocità sulla ventola 1 segnala un arresto inaspettato.	Spegnere e riavviare il dispositivo. Se il messaggio d'errore persiste, chiamare un tecnico del servizio di assistenza.	•
8	La misurazione della velocità sulla ventola 2 segnala un arresto inaspettato.	Spegnere e riavviare il dispositivo. Se il messaggio d'errore persiste, chiamare un tecnico del servizio di assistenza.	•
9	La misurazione della velocità sulla ventola 3 segnala un arresto inaspettato.	Spegnere e riavviare il dispositivo. Se il messaggio d'errore persiste, chiamare un tecnico del servizio di assistenza.	•2
10	La misurazione della velocità sulla ventola 4 segnala un arresto inaspettato.	Spegnere e riavviare il dispositivo. Se il messaggio d'errore persiste, chiamare un tecnico del servizio di assistenza.	•2
12	Errore di avvio azionamento - segnale di velocità non ricevuto.	La piattaforma dell'agitatore è bloccata. Controllare se lo spazio intorno alla piattaforma è sufficiente e/o ridurre il carico (accessori e campioni) sulla piattaforma. Successivamente, premere un'altra volta START sull'agitatore. Se il messaggio d'errore persiste, contattare un tecnico del servizio di assistenza.	٩
24	Errore di misurazione velocità durante un ciclo.	È stato rilevato un cambiamento anomalo della velocità. Controllare il carico (accessori e campioni) e/o i supporti sulla piattaforma dell'agitatore. Successivamente, premere un'altra volta START sull'agitatore. Se il messaggio d'errore persiste, contattare un tecnico del servizio di assistenza.	
26	Misurazione velocità ha rilevato un fermo inaspettato durante un ciclo.	Assicurarsi che lo spazio intorno alla piattaforma sia sufficiente e/o ridurre il carico (accessori e campioni) sulla piattaforma. Assicurarsi che il fusibile (interruttore automatico) sul retro dell'agitatore non sia scattato («Fusibile» a pagina 48). Successivamente, premere un'altra volta START sull'agitatore. Se il messaggio d'errore persiste, contattare un tecnico del servizio di assistenza.	٢
82	La misurazione della corrente motore ha rilevato un sovraccarico.	È stata rilevata una sovracorrente motore. Non caricare o scaricare la piattaforma durante il funzionamento. Ridurre la velocità o adattare il carico (accessori e campioni) sulla piattaforma.	— 9
83	Misurazione della corrente motore fuori dai limiti.	È stata rilevata una sovracorrente motore. Non caricare o scaricare la piattaforma durante il funzionamento. Ridurre la velocità o adattare il carico (accessori e campioni) sulla piattaforma.	— 0

Tabella 59: Lista degli errori, avvertimenti e allarmi

AVVISO Se viene visualizzato un messaggio d'errore non contenuto in questa tabella, contattare un tecnico del servizio di assistenza.

GPL (Licenza Pubblica Generica)

Parti del software del dispositivo usano un software open-source, pubblicato sotto la licenza GPL, LGPL o un'altra licenza open-source. In dettaglio, si tratta delle librerie elencate nella tabella. Il codice sorgente delle librerie usate (librerie di terzi) può essere ottenuto da Thermo Fisher Scientific se la rispettiva licenza lo prevede. Le rispettive condizioni di licenza del software open-source usato fanno parte del pacchetto di codice sorgente messo a disposizione.

Librerie di terzi

Libreria	Versione	Impiego via	Licenza
Qt	5.8	BSP	LGPLv3
Log4Cplus	1.2.0	Applicazione	Apache Public License v2 / Licenza BSD con due clausole
boost	1.72.0	Applicazione	Licenza Boost 1.0
json11	1.0.0	Applicazione	Licenza MIT

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

Indice analitico

Α

Accessori 18, 89 Allarmi 70 Autoclavaggio 126 Avvertimenti 70

С

Calibrazione 73 Calibrazione di temperatura 127 Caricamento 116, 117 Collegamento alla rete 48 Componenti di metallo 124 Componenti in plastica 124 Configurazione iniziale 55 Controllo dell'accesso 72 Cura 122

D

Data 75 Dati tecnici 12 Decontaminazione 125 Diagrammi 87 Direttive 44 Disinfezione 125 Display 74 Dotazione di fornitura 50 Durata della tacitazione temporanea (Snooze) 71

Ε

Elementi di comando 74 Eliminazione guasti 136 Ethernet 48

F

File & Info 78 Formato data 76 Funzionamento 61 Fusibile 48

G

GPL (Licenza Pubblica Generica) 138

Impostazione della temperatura 65 Impostazioni 69 Informazioni generali sul prodotto 45 Installazione dei rack portaprovette 97 Installazione dei supporti ad angolo regolabile per rack portaprovette 98 Installazione dei supporti per micropiastra / piastra DeepWell 99 Installazione dei supporti porta matracci e dei recipienti 93 Installazione dei supporti quadrati 95 Installazione dei supporti regolabili per recipienti 100 Installazione del collettore del gas 114 Installazione della piattaforma 90 Installazione del rack porta becher 102 Installazione del supporto multiuso 106 Installazione del supporto per imbuto separatore 103 Installazione e uso del tappetino antiscivolo 107 Interfaccia utente grafica 61 Ispezione degli accessori 124

L

Lingua 75 Luminosità 74 Luogo di installazione 51

Μ

Manutenzione 78, 122, 135 Modalità tempo 74 Modo veglia 76

Ν

Nome dell'unità 77 Norme 44

Ρ

Parole e simboli di segnalazione 6 Personalizzazione del menu 77 Piattaforme 19 Piattaforme a due livelli 92 Piattaforme universali 91 Programmi 79 Pulizia 124

R

Raccordi 47 Regione 77 Registro degli eventi 85 Riavvio automatico 74

S

Simboli utilizzati nelle istruzioni 7 Smaltimento 135 Specifiche tecniche 12 Spedizione 60, 135

Т

Temp nominale 65, 66 Tempo 76 Trasporto 53

U

USB 49 Uso previsto 5, 116, 118

V

Valore nominale della temperatura 74 Valore nominale di velocità 74 Velocità massima 119 Velocità sicura 118 Volume 70

thermo scientific



Prodotto per

Thermo Electron LED GmbH Zweigniederlassung Osterode Am Kalkberg, 37520 Osterode am Harz Germany Nazione di origine: 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 è il manuale d'uso originale. Questo manuale di istruzioni è una traduzione del manuale di istruzioni originale.

thermofisher.com/shaker

© 2019–2024 Thermo Fisher Scientific Inc. Tutti i diritti riservati.

Se non specificato diversamente, tutti i marchi di fabbrica sono proprietà della Thermo Fisher Scientific Inc. e delle sue controllate. Non tutti i prodotti sono disponibili in tutti i paesi. Per maggiori dettagli consultare il rappresentante commerciale locale.

Le immagini pubblicate all'interno delle presenti istruzioni per l'uso servono unicamente come riferimento. Le impostazioni e le lingue illustrate possono differire.

Australia +61 39757 4300

Austria +43 1 801 40 0

Belgio +32 53 73 42 41

Cina +800 810 5118 **o** +400 650 5118

Francia +33 2 2803 2180

Germania nazionale, numero verde 0800 1 536 376

Germania internazionale +49 6184 90 6000 India +91 22 6716 2200

Italia +39 02 95059 552

Giappone +81 3 5826 1616

Paesi Bassi +31 76 579 55 55

Nuova Zelanda +64 9 980 6700

Paesi nordici/del Baltico/CSI +358 10 329 2200

Russia +7 812 703 42 15 **Spagna/Portogallo** +34 93 223 09 18

Svizzera +41 44 454 12 12

Gran Bretagna/Irlanda +44 870 609 9203

Stati Uniti / Canada +1 866 984 3766

Altri paesi asiatici +852 2885 4613

Altri paesi +49 6184 90 6000

