

Thermo Scientific Solarisシェーカー(振とう機) シリーズ

取扱説明書

70900201-е • 2024-11



目次

前書き	5
取扱説明書概要	5
定められた用途..........................	5
開放型シェーカー	5
恒温シェーカー(加熱式)および冷却シェーカー	
(冷却・加熱式)	5
操作適格者	5
注意喚起語とシンボル.....................	6
シェーカーと付属品に使用されるシンボル	6
取扱説明書で使用されるシンボル	7
安全上の注意...........................	7
1 技術仕様 1	1
	1
$I. I. \text{Therma Scientific Scleric } (\lambda = 1.7) 2000$	I
・ 「	1
Thermo Scientific Solaris (ソラリス)	
4000 シェーカー(振とう機)	2
Thermo Scientific Solaris 2000 シェーカー1	3
Thermo Scientific Solaris 4000 I シェーカー1	4
Thermo Scientific Solaris 2000 R シェーカー1	5
Thermo Scientific Solaris 4000 R シェーカー1	6
1. 2. 付属品	7
1. 2. 1. プラットフォーム(振とう台) 1	8
1. 2. 2. フラスコクランプ 1	9
1. 2. 3. マイクロプレート・ディープウェルプレートクランプ 2	23
1. 2. 4. 試験管立て	:4
1. 2. 5. 角度調節式試験管立てホルダー	8
1. 2. 6. 調節式容器クランプ	0
1. 2. 7. 角型メディアクランプ 3	2
1. 2. 8. Nalgeneビーカーラック 3	4
1. 2. 9. 分液ロートクランプ 3	8
1. 2. 10. ユーティリティトレイ	9

	1. 2	. 11.	粘着マット	40
	1. 2	. 12.	その他の付属品...................	42
1.	3.	指令	および 規格	43
1.	4.	製品	概要...................	44
	1. 4	. 1.	Solaris 2000	44
	1. 4	. 2.	Solaris 4000	44
	1. 4	. 3.	Solaris 2000 I / 2000 R	45
	1. 4	. 4.	Solaris 4000 I / 4000 R	45
	1. 4	. 5.	接続	46
2.	運搬	と設調	置	49
2.	1.	開梱		49
		納	品内容	49
2.	2.	設置	場所	50
2	3.	運搬		52
	•	~シ	· エーカーの取扱い	52
2.	4.	水平	調整方法.....................	53
2.	5.	電源		53
2	6.	初回	記動	54
2	7	保管		59
2.	γ. Q	() () () () () () () () () () () () () (50
	0.	廷公		
3.	運転			60
3.	1.	電源	のオン/オフ	60
3.	2.	グラ	フ表示のあるユーザーインターフェース	60
	3. 2	. 1.	シェーカーの基本操作パラメーターの設定....	61
	3. 2	. 2.	ステータス	66
	3. 2	. 3.	セッティング	68
	3. 2	. 4.	プログラム.................	78
3.	3.	付属	品	88
	3. 3	. 1.	プラットフォームの設置............	89
	3. 3	. 2.	フラスコクランプと容器の取付け.......	92
	3. 3	. 3.	角型クランプの取付け.............	94
	3. 3	. 4.	試験管立ての取付け..............	96
	3. 3	. 5.	角度調節式試験管立てホルダーの取付け.....	97

	3. 3. 6. マイクロプレート・ディープウェルプレートの クランプの取付け	00
	クランノの取り1、	90
	3 3 8 ビーカーラックの取付け	101
	3. 3. 9. 分液ロートクランプの取付け	102
	3. 3. 10. ユーティリティトレイの取付け	105
	3. 3. 11. 粘着マットの取付けと使用	106
	3. 3. 12. ガス供給マニホールドの取付け	113
3.	4. 負荷と 通常操作	115
	読み込む......................	116
	通常操作	117
4.	メンテナンスとケア................	121
4.	1. 基本事項	122
	付属品の点検......................	123
4.	2. クリーニング	123
	タッチスクリーン	123
4.	3. 消毒	124
4.	4. 除染	124
4.	5. オートクレーブ	125
4.	6. 温度校正	126
	4. 6. 1. 温度校正の最良実践法	126
	4. 6. 2. 温度の校正の手順	126
	4. 6. 3. 温度校正の探索	130
4.	7. ファームウェアのインストール	130
4.	8. プラットフォームの交換	134
4.	9. アフターサービス	134
4.	10. 送付と廃棄	134
5.	トラブルシューティング..............	135
GPL	(General Public License:一般公有使用許諾)	137

前書き

取扱説明書概要

本取扱説明書では、下記のシェーカーに対応している付属品 (17ページの 「1.2. 付属品」)について説明しています。

シェーカー	製品番号
Thermo Scientific Solaris 2000 $\succ \pm - \pi -,$ 100–240 V ± 10 %, 50 / 60 Hz	SK2000
Thermo Scientific Solaris 4000 \not x — \not — $,$ 100–240 V ±10 %, 50 / 60 Hz	SK4000
Thermo Scientific Solaris 2000 I $\mathrel{\vartriangleright} \mathtt{x}-\mathtt{h}-$, 100–120, 200–240 V ± 10 %, 50 / 60 Hz	SK2001
Thermo Scientific Solaris 4000 I $\mathrel{\vartriangleright} \pm \pm - \pm -$, 100–120, 200–240 V ± 10 %, 50 / 60 Hz	SK4001
Thermo Scientific Solaris 2000 R $\mathrel{\vartriangleright} \pm \pm - \pm -$, 100–240 V ± 10 %, 50 / 60 Hz	SK2002
Thermo Scientific Solaris 4000 R $\succ\pm-\pi-$, 100-240 V ±10 %, 50 / 60 Hz	SK4002

本書に記載されている説明と手順には、製品名を明記する代わりに、以下のよう な分類別にシェーカーが記載されています。

- 開放型シェーカー
- 恒温シェーカー(加熱式)
- 冷却シェーカー(冷却・加熱式)

定められた用途

開放型シェーカー

アプリケーションを用いて速度と時間を制御・管理のもと、2次元軌道で溶液を 振とうさせるための、研究および製造目的に使用される装置です。

恒温シェーカー(加熱式)および冷却シェーカー(冷却・加熱式) アプリケーションを用いて速度、時間、温度を制御・管理のもと、2次元軌道で 溶液を振とうさせるための、研究および製造目的に使用される装置です。

操作適格者

このシェーカーのご利用は操作訓練を受けた人に限ります。

注意喚起語とシンボル

注意喚起語	危険の度合い
警告	これを避けなければ、死亡や重症の怪我を負う可能性のある、危険 な状況を示します。
注意	これを避けなければ、中軽傷の怪我を負う可能性のある、危険な状 況を示します。
注	重要な情報を示していますが、危険に関する物ではありません。

シェーカーと付属品に使用されるシンボル

安全にお使いいただくために必ず取扱説明書の記載事項に従ってください。

8	取扱説明書参照
	電源プラグを抜く
	一般的な危険
	高温表面の接触による危険性
	機械部品が閉まる事故によって指 や手が負傷する危険性

取扱説明書で使用されるシンボル

ご使用の際は、周囲を危険にさらさぬよう、取扱い上の指示をお守りください。

一般的な危険	<u>A</u>	電気的危険性
生物学的危険性		切り傷を負う危険性
可燃物による危険性	i	重要な情報を示していま すが、危険に関する物で はありません。
機械部品が閉まる事故に よって指や手が負傷する 危険性。		

安全上の注意



警告

これらの安全上の注意に従わない誤使用は死亡または重傷を 伴なう危険性を誘発する場合がを誘発する場合があります。

- 安全上の注意をお守りください。
- シェーカー本来の用途以外では使用しないでください。
 誤った使用方法は破損、汚染、および生命に関わる怪我の原因となる恐れがあります。
- このシェーカーのご利用は操作訓練を受けた人に限ります。
- オペレーターには適切な個人用保護具使用の確認義務が あります。世界保健機関(WHO)の実験室バイオセーフ ティ指針、ならびに使用国における規定にご注意ください。



不適切な電源使用による破損。 シェーカーが適切に接地されたソケットに差し込まれている ことを確認してください。

警告



警告

有害物質の取り扱いによる危険。 塩類溶液、酸、塩基などの腐食性のサンプルを取り扱ったと きは、必ず付属品とシェーカーを十分に清掃してください。

- シェーカーは不活性機器でも防爆機器でもありません。
 発火しやすい場所では絶対に使用しないでください。
- 有毒物質や放射性物質、あるいは病原微生物類の振とう
 には、適切な安全対策を施してください。
- 有害物質を振とうさせる際は、世界保健機関(WHO)の 「実験室バイオセーフティマニュアル」および現地の規 定を遵守してください。世界保健機関(WHO)の「実験 室バイオセーフティマニュアル」に従って、リスク群2 の微生物サンプルを振とうさせる場合は、エアロゾル封 じ込め用バイオロジカルシールを必ず使用して用バイオ ロジカルシールを必ず使用してください。世界保健機関 WHOのホームページ (www.who.int) にて、「実験室バイ オセーフティ指針」をご検索ください。価の高いリスク グループの物質の場合、予防対策を必ず二つ以上備えて ご利用ください。
- シェーカーやシェーカーの部品が毒素や病原物質に汚染 された場合は、適切な消毒および除染対策を取る必要が あります(124ページの「除染」、124ページの「消 毒」の参照)。
- 危険な状況が発生した場合は、シェーカーの電源を切り、速やかにその場を離れてください。



警告

感染性物質による健康被害。

事故による流出などで液体や物がプラットフォームの下に入り込んだ場合は、直ちにシェーカーの電源を切り、プラグを 抜き、プラットフォームを取り外してください(18ページの「プラットフォーム(振とう台)」)。

通常の検査法に従って、流出物を清掃してください。適切な 個人用保護具を使用してください。



爆発性または可燃性の物質や素材の振とうによる健康被害。 爆発性や可燃性のある物質や素材はシェーカーにかけないで ください。

警告



高温表面の接触による火傷。

注意

温度調節機能付シェーカーでは、フードの内側で温度が上昇 し、プラットフォームや付属品などの表面が熱くなる可能性 があります。高温になっている表面に触れないでください。 表面が高温になっている場合は冷えるまで待ちます。



注意

ガラス破片による切り傷。

プラットフォームやクランプなどの付属品が外れると、容器 がシェーカーから落ちてガラスが割れることがあります。

プラットフォームと付属品が正しい工具とネジで適切に取り 付けられているか確認してください。クランプが、選択され た速度で容器とサンプルの負荷に耐えることができるか確認 してください。個別に設定された負荷のクランプがシェーカ ーの安定限界速度よりも遅いことがあります。115ページの 「3.4.負荷と 通常操作」してください。

異常な音に注意してください。異常な音が聞こえる場合は、 プラットフォームや付属品が外れている可能性があります。



注意

- タッチスクリーン破損によるシェーカーの損傷・誤動作。
- タッチスクリーンが破損している場合は、シェーカーを 操作しないでください。
- シェーカーの電源を切り、電源プラグを抜いてください。認定を受けたサービス技術者にタッチスクリーンの 交換を要請してください。



注意

容器の破損・液漏れによる生物学的危害。

付属品が正しく取り付けられていないとサンプルが流出する 可能性があります。

- 付属品が正しい工具とネジで適切に取り付けられている か確認してください。
- 付属品がプラットフォームに適切に組み合わさっている か確認してください。
- 容器は必ず適切な寸法の付属品で使用してください。
- 容器は損傷がないものを正しく取り付けてください。



装填の不備や付属品の損傷は安全性を低下させる恐れがあり ます。

注意

- 必ず負荷(付属品とサンプル)ができるだけ均等に分散 されるように設置してください。特にデュアルスタック プラットフォームを操作する場合は注意してください。
- 腐食や亀裂の兆しが認められる付属品は使用しないでく ださい。詳細については、カスタマーサービスにお問い 合わせください。
- 適切に装填されたシェーカーのみ使用してください。
- シェーカーに過大な負荷をかけないでください。
- シェーカーを操作する前に、付属品が正しく取り付けられているか確認してください。88ページの「付属品」の指示に従ってください。



注意

作業基礎の無視に起因する身体的危害。

- プラットフォームが適切に設置されていない状態でシェ
 ーカーを操作しないでください。
- 部品の外装が損傷または欠損している場合は、シェーカ ーを使用しないでください。
 - フードのスプリングに不具合のあるシェーカーは使用しないでください。フードのスプリングに不具合がある場合は、いっぱいまで開けた位置でフードを安全に保持することができません。
 - シェーカーの作動中は、シェーカーを移動させないでく ださい。
- シェーカーに寄り掛かからないでください。
- シェーカーが完全に停止し、タッチスクリーンで停止したことを確認するまではシェーカーに物を載せたり降ろしたりしないでください。
- シェーカーの作動中は上に物を置かないでください。
- シェーカーの作動中はプラットフォームや付属品に触れ ないでください。
- オペレーターはシェーカーのハウジングを開けてはいけません。
- 適合しない付属品の使用は損傷防止機能を損なうことがあり ます。

注

1

このシェーカーにはThermo Fisher Scientificが承認し た付属品のみご利用ください。最新の製品リストについて は、www.thermofisher.comをご覧ください。



注

シェーカーを停止するには以下の操作を行います。 STOPキーを押します。シェーカーのメインスイッチで電源を 切ります。電源コンセントを引き抜きます。緊急時には、電 源を切断してください。

10 / 138

1. 技術仕様

1. 1. 仕様

Thermo Scientific Solar シェーカー(振とう機)	ris (ソラリス	K) 2000	
速度範囲	15~525rpm		
実行時間	99時間59分(1分刻み)	または連続モード	
最高速度時の騒音レベル	50dB(A)(装置正面加	vら1 m、高さ1.6m)	
最大負荷重量(プラットフォーム、付属 品、サンプルを含む)	25kg		
電源	100-240 V ±10 %, 50	/ 60 Hz	
消費電力	60 W		
環境条件			
保管・出荷時	温度: -10 °C ~ 55 湿度: 15%~85%	°C	
操作時	屋内使用 標高海抜3000 mまで 温度:5°C~40°C 31°C以下で最高相対語 に減少	星度80%40°Cで相対湿度	50%まで一次関数的
汚染度	2		
過電圧カテゴリ	II		
IPコード (保護等級)	20		
インターフェース	USB	イーサネット	
	2x USB-A 2.0	RJ45	
寸法	長さ	中国	高さ
	47 cm	37 cm	15 cm
重量	20. 9kg		

表 1: 技術データ Solaris 2000

Thermo Scientific Solaris (ソラリス) 4000 シェーカー (振とう機)					
速度範囲	15~525rpm				
実行時間	99時間59分(1分刻み)) または連続モード			
最高速度時の騒音レベル	54dB(A)(装置正面)	から1 m、高さ1.6m)			
最大負荷重量(プラットフォーム、付属 品、サンプルを含む)	43kg				
電源	100-240 V ±10 %, 50	0 / 60 Hz			
消費電力	80 W				
環境条件					
保管・出荷時	温度: −10 °C ~ 55 湿度: 15%~85%	°C			
操作時	屋内使用 標高海抜3000 mまで 温度: 5°C~40°C 31°C以下で最高相対; に減少	显度80%40°Cで相対湿度	50%まで一次関数的		
汚染度	2				
過電圧カテゴリ	II				
IPコード(保護等級)	20				
インターフェース	USB	イーサネット			
	2x USB-A 2.0	RJ45			
寸法	長さ	中畐	高さ		
	65 cm	58 cm	18 cm		
重量	75. 1kg				

表 2: 技術データ Solaris 4000

Thermo Scientific Sola	ris 2000 シュ	∟ −カー	
速度範囲	15~525rpm		67.
温度設定幅	30-60 ° C		
フラスコ内の温度安定性 37 °Cの場合	周囲温度23°C、安定し ± 0.1 °C	ンたチャンバー温度で1時	寺間
フラスコ内の温度均一性 37 °Cの場合	周囲温度23°C、安定し ± 0.5 °C	、たチャンバー温度で1時	节 目
実行時間	99時間59分(1分刻み)	または連続モード	
最高速度時の騒音レベル	52dB(A)(装置正面カ	ヽら1 m、高さ1.6m)	
最大負荷重量(プラットフォーム、付属 品、サンプルを含む)	16kg (35 lb) 2.3kg (5.1 lb) のプラ を含む	ラットフォーム (11x14a	インチ : 約28x36cm)
電源	100-120, 200-240 V ±	=10 %, 50 / 60 Hz	
消費電力	900 W		
環境条件			
保管・出荷時	温度: -10 °C ~ 55 ′ 湿度: 15%~85%	°C	
操作時	屋内使用 標高海抜3000 mまで 温度: 5°C〜40°C 31°C以下で最高相対造 に減少	2度80%40°Cで相対湿度	50%まで一次関数的
汚染度	2		
過電圧カテゴリ	II		
IPコード(保護等級)	20		
インターフェース	USB	イーサネット	
	2x USB-A 2.0	RJ45	
寸法	長さ	幅	高さ
	70 cm	36 cm	46 cm
フードを開けた状態			79 cm
重量	44. 9kg		

表 3: 技術データ Solaris 2000 I

Thermo Scientific Solar	is 4000 I S	ンェーカー			
速度範囲	15~525rpm				
温度設定幅 30-60 ° C					
フラスコ内の温度安定性 37 ° Cの場合	周囲温度23°C、安定し ± 0.1 °C	ったチャンバー温度で1時	[]]		
フラスコ内の温度均一性 37 ° Cの場合	周囲温度23°C、安定し ± 0.5 °C	ったチャンバー温度で1時	- 月月 月 月		
実行時間	99時間59分(1分刻み)	または連続モード			
最高速度時の騒音レベル	52dB(A)(装置正面か	vら1 m、高さ1.6m)			
最大負荷重量(プラットフォーム、付属 品、サンプルを含む)	32kg (70.51b) 4.3kg (9.51b) のプラ を含む	ットフォーム(18x18イ)	ンチ : 約46x46cm)		
電源	100-120, 200-240 V ±	±10 %, 50 / 60 Hz			
消費電力	900 W				
環境条件					
保管・出荷時	温度: -10 °C ~ 55 湿度: 15%~85%	°C			
操作時	屋内使用 標高海抜3000 mまで 温度:5°C~40°C 31°C以下で最高相対語 に減少	星度80%40°Cで相対湿度8	50%まで一次関数的		
汚染度	2				
過電圧カテゴリ	II				
IPコード(保護等級)	20				
インターフェース	USB	イーサネット			
	2x USB-A 2.0	RJ45			
寸法	長さ	中国	高さ		
	77 cm	57 cm	55 cm		
フードを開けた状態			94 cm		
重量	68. 1kg				

表 4: 技術データ Solaris 4000 I

Thermo Scientific Sola	ris 2000 R 🗄	シェーカー	
速度範囲	15~525rpm		C .
温度設定幅	5-60 ° C		
フラスコ内の温度安定性 37 °Cの場合	周囲温度23°C、安定し ± 0.1 °C	ったチャンバー温度で1時	寺間
フラスコ内の温度均一性 37 °Cの場合	周囲温度23°C、安定し ± 0.5 °C	したチャンバー温度で1時	寺間
実行時間	99時間59分(1分刻み)	または連続モード	
最高速度時の騒音レベル	56dB(A)(装置正面 <i>f</i>	から1 m、高さ1.6m)	
最大負荷重量(プラットフォーム、付属 品、サンプルを含む)	16kg (35 lb) 2.3kg (5.1 lb) のプラ を含む	ラットフォーム(11x14a	インチ : 約28x36cm)
電源	100-240 V ±10 %, 50	/ 60 Hz	
消費電力	350 W		
環境条件			
保管・出荷時	温度: -10 °C ~ 55 湿度: 15%~85%	°C	
操作時	屋内使用 標高海抜3000 mまで 温度: 5°C~40°C 31°C以下で最高相対語 に減少	显度80%40°Cで相対湿度	50%まで一次関数的
汚染度	2		
過電圧カテゴリ	II		
IPコード(保護等級)	20		
インターフェース	USB	イーサネット	
	2x USB-A 2.0	RJ45	
寸法	長さ	中国	高さ
	70 cm	36 cm	46 cm
フードを開けた状態			79 cm
重量	47. 2kg		

表 5: 技術データ Solaris 2000 R

Thermo Scientific Sola	ris 4000 R ジ	ィェーカー			
速度範囲	15~525rpm				
温度設定幅	4-60 ° C				
フラスコ内の温度安定性 37 °Cの場合	周囲温度23°C、安定し ± 0.1 °C	たチャンバー温度で1時	間		
フラスコ内の温度均一性 37 °Cの場合	周囲温度23°C、安定し ± 0.5 °C	周囲温度23°C、安定したチャンバー温度で1時間 ± 0.5 °C			
実行時間	99時間59分(1分刻み)	または連続モード			
最高速度時の騒音レベル	56dB (A) (装置正面か	ら1 m、高さ1.6m)			
最大負荷重量(プラットフォーム、付属 品、サンプルを含む)	32kg (70.51b) 4.3kg (9.51b) のプラ、 を含む	ットフォーム(18x18イン	ンチ:約46x46cm)		
電源	100-240 V ±10 %, 50	/ 60 Hz			
消費電力	750 W				
環境条件					
保管・出荷時	温度: -10 °C ~ 55 ° 湿度: 15%~85%	С			
操作時	屋内使用 標高海抜3000 mまで 温度: 5°C~40°C 31°C以下で最高相対湿 に減少	度80%40° Cで相対湿度5	0%まで一次関数的		
汚染度	2				
過電圧カテゴリ	II				
IPコード(保護等級)	20				
インターフェース	USB	イーサネット			
	2x USB-A 2.0	RJ45			
寸法	長さ	中国	高さ		
	77 cm	57 cm	55 cm		
フードを開けた状態			94 cm		
重量	74. 9kg				

表 6: 技術データ Solaris 4000 R

1. 2. 付属品



適合しない付属品の使用は損傷防止機能を損なうことがありま す。

このシェーカーにはThermo Fisher Scientificが承認した付属 注意 品のみご利用ください。

最新の製品リストについては、www.thermofisher.comをご覧ください。

1. 2. 1. プラットフォーム(振とう台)

プラットフォーム		
Solaris 2000	重量	7
Thermo Scientific Solaris 12x14 標準ブラットフォーム	2. 7kg	SK1214
Thermo Scientific Solaris 12x14 デュアルスタック標準プラットフォーム	6. 4kg	SK1214D
Thermo Scientific Solaris 12x14 デュアルスタック標準プラットフォ ーム・アップグレード用キット		SK1214DK
Thermo Scientific Solaris 18x18 標準ブラットフォーム	4. 3kg	SK1818
Thermo Scientific Solaris 18x18 デュアルスタック標準プラットフォーム	9.7kg	SK1818D
Thermo Scientific Solaris 18x18 デュアルスタック標準プラットフォ ーム・アップグレード用キット		SK1818DK
Thermo Scientific Solaris 18x24 標準ブラットフォーム	6. 0kg	SK1824
Solaris 4000		
Thermo Scientific Solaris 18x30 標準ブラットフォーム	7. 5kg	SK1830
Thermo Scientific Solaris 18x30 デュアルスタック標準プラットフォーム	16. 3kg	SK1830D
Thermo Scientific Solaris 18x30 デュアルスタック標準プラットフォ ーム・アップグレード用キット		SK1830DK
Thermo Scientific Solaris 36x24 標準プラットフォーム	12. 0kg	SK3624
Solaris 2000 I / 2000 R		
Thermo Scientific Solaris 11x14 標準プラットフォーム	2. 3kg	SK1114
Solaris 4000 I / 4000 R		
Thermo Scientific Solaris 18x18 標準ブラットフォーム	4. 3kg	SK1818
スペアキットと付属品		
クランプスペアキット (ネジ)		SK1001
プラットフォームスペアキット、小(プラットフォーム用ネジ、工具)		SK0100
ブラットフォームスペアキット、大(プラットフォーム用ネジ、工具)		SK0101
付属品用ドライバー		75004131

表 7: プラットフォーム製品

1. 2. 2. フラスコクランプ



<u>Solaris 2000プラットフォーム用</u>

	製品番号	標準		デュアルスタック		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
エルレンマイヤー フラス コ 10ml	30150BI	72	113	157	140	226
エルレンマイヤー フラス コ 25ml	30151	42	64	80	80	124
エルレンマイヤー フラス コ 50ml	30152BI	42	64	80	80	124
エルレンマイヤー フラス コ 125ml	30153	15	32	40	30	62
エルレンマイヤー フラス コ 250ml	30154BI	9	16	24	16	32
エルレンマイヤー フラス コ 300ml	30155	9	16	20	16	32
エルレンマイヤー フラス コ 500ml	30156BI	9	16	20	16	32
エルレンマイヤー フラ スコ 11	30157BI	4	9	11	8	16
エルレンマイヤー フラ スコ 21	30158	3	5	6	-	-
エルレンマイヤー フラ スコ 41	30159	1	4	4	-	-
エルレンマイヤー フラ スコ 51	30159B	1	2	4	-	-
エルレンマイヤー フラ スコ 61	30160	1	2	2	-	-
フェルンバッハ フラスコ 2,800ml	30162	1	4	4	-	-
Low-Form培養フラスコ 2.5	30161	1	2	2	-	-

表 8: Solaris 2000プラットフォーム用フラスコクランプ製品

<u>Solaris 4000プラットフォーム用</u>

	製品番号	Ŕ	票準	デュアルスタ ック
		18 x 30	36 x 24	18 x 30
エルレンマイヤー フラス コ 10ml	30150BI	203	187	402
エルレンマイヤー フラス コ 25ml	30151	112	187	220
エルレンマイヤー フラス コ 50ml	30152BI	112	187	220
エルレンマイヤー フラス コ 125ml	30153	46	83	92
エルレンマイヤー フラス コ 250ml	30154BI	34	40	68
エルレンマイヤー フラス コ 300ml	30155	28	40	56
エルレンマイヤー フラス コ 500ml	30156BI	28	40	56
エルレンマイヤー フラ スコ 11	30157BI	14	20	28
エルレンマイヤー フラ スコ 21	30158	8	11	16
エルレンマイヤー フラ スコ 41	30159	6	8	-
エルレンマイヤー フラ スコ 51	30159B	5	8	-
エルレンマイヤー フラ スコ 61	30160	3	6	-
フェルンバッハ フラスコ 2,800ml	30162	6	8	12
Low-Form培養フラスコ 2.5	30161	3	6	6

表 9: Solaris 4000プラットフォーム用フラスコクランプ製品

	製品番号	標準
		11 x 14
エルレンマイヤー フラスコ 10ml	30150BI	59
エルレンマイヤー フラスコ 25ml	30151	35
エルレンマイヤー フラスコ 50ml	30152BI	35
エルレンマイヤー フラスコ 125ml	30153	15
エルレンマイヤー フラスコ 250ml	30154BI	7
エルレンマイヤー フラスコ 300ml	30155	7
エルレンマイヤー フラスコ 500ml	30156BI	7
エルレンマイヤー フラスコ 11	30157BI	4
エルレンマイヤー フラスコ 21	30158	2
エルレンマイヤー フラスコ 41	30159	-
エルレンマイヤー フラスコ 51	30159B	-
エルレンマイヤー フラスコ 61	30160	-
フェルンバッハ フラスコ 2,800ml	30162	1
Low-Form培養フラスコ 2.5	30161	1

<u>Solaris 2000 I / 2000 Rプラットフォーム用</u>

表 10: Solaris 2000 I/2000 Rプラットフォーム用フラスコクランプ製品

	製品番号	標準
		18 x 18
エルレンマイヤー フラスコ 10ml	30150BI	113
エルレンマイヤー フラスコ 25ml	30151	64
エルレンマイヤー フラスコ 50ml	30152BI	64
エルレンマイヤー フラスコ 125ml	30153	32
エルレンマイヤー フラスコ 250ml	30154BI	16
エルレンマイヤー フラスコ 300ml	30155	16
エルレンマイヤー フラスコ 500ml	30156BI	16
エルレンマイヤー フラスコ 11	30157BI	9
エルレンマイヤー フラスコ 21	30158	6
エルレンマイヤー フラスコ 41	30159	-
エルレンマイヤー フラスコ 51	30159B	-
エルレンマイヤー フラスコ 61	30160	-
フェルンバッハ フラスコ 2,800ml	30162	4
Low-Form培養フラスコ 2.5	30161	1

<u>Solaris 4000 I / 4000 Rプラットフォーム用</u>

表 11: Solaris 4000 I/4000 Rプラットフォーム用フラスコクランプ製品

2. 3. マイクロプレート・
 ディープウェルプレートクランプ



<u>Solaris 2000プラットフォーム用</u>

	製品番 号		標準		デュアル	スタック
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
マイクロプレート・ディー プウェルプレート	30175	7	12	16	14	24

表 12: Solaris 2000プラットフォーム用マイクロプレート・ディープウェルプレートクラ ンプの容量

<u>Solaris 2000 I / 2000 Rプラットフォーム用</u>

	製品番号	標準
		11 x 14
マイクロプレート・ディー プウェルプレート	30175	6

表 13: Solaris 2000 I/2000 Rプラットフォーム用マイクロプレート・ディープウェルプ レートホルダーの容量

<u>Solaris 4000プラットフォーム用</u>

	製品番号	標準		デュアルスタ ック
		18 x 30	36 x 24	18 x 30
マイクロプレート・ディー プウェルプレート	30175	23	36	42

表 14: Solaris 4000プラットフォーム用マイクロプレート・ディープウェルプレートホル ダーの容量

<u>Solaris 4000 I / 4000 Rプラットフォーム用</u>

	製品番号	標準
		18 x 18
マイクロプレート・ディー プウェルプレート	30175	10

表 15: Solaris 4000 I/4000 Rプラットフォーム用マイクロプレート・ディープウェルプ レートホルダーの容量

23 / 138

1. 2. 4. 試験管立て

<u>Solaris 2000プラットフォーム用</u>



	製品番号		標準		デュアル	スタック
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
ハーフサイズ						
10~13mm, 赤、配列 6x6	30181	8	12	15	14	22
14~16mm, オレンジ、配列 6x6	30183	5	9	11	9	16
17~20mm, 白、配列 4x5	30185	7	11	14	12	20
21~25mm、 青、配列 4x4	30187	6	9	11	10	16
26~30mm、 緑、配列 3x3	30189	6	9	12	11	17
1.5ml 微量遠心機用、 青、配列 4x6	30191	6	10	13	11	19
フルサイズ						
10~13mm、配列 6x12	30180BI	3	7	9	6	12
14~16mm、配列 6x12	30182	3	4	6	6	8
17~20mm、配列 4x10	30184	3	5	7	6	8
21~25mm、配列 4x10	30186	2	3	5	4	6
26~30mm、配列 3x8	30188	3	4	5	5	8
1.5ml 微量遠心機用、 配列 8x12	30190	3	4	7	6	8

表 16: Solaris 2000プラットフォーム用試験管立て製品

<u>Solaris 4000プラットフォーム用</u>

	製品番号	桴	票準	デュアルスタック
		18 x 30	36 x 24	18 x 30
ハーフサイズ				
10~13mm, 赤、配列 6x6	30181	21	32	40
14~16mm, オレンジ、配列 6x6	30183	13	20	24
17~20mm, 白、配列 4x5	30185	18	20	34
21~25mm、 青、配列 4x4	30187	14	22	26
26~30mm、 緑、配列 3x3	30189	15	24	30
1.5ml 微量遠心機用、 青、配列 4x6	30191	18	24	34
フルサイズ				
10~13mm、配列 6x12	30180BI	13	20	18
14~16mm、配列 6x12	30182	9	12	16
17~20mm、配列 4x10	30184	9	15	18
21~25mm、配列 4x10	30186	6	9	12
26~30mm、配列 3x8	30188	7	10	13
1.5ml 微量遠心機用、 配列 8x12	30190	9	12	17

表 17: Solaris 4000プラットフォーム用試験管立て製品

25 / 138

	製品番号	標準
		11 x 14
ハーフサイズ		
10~13mm, 赤、配列 6x6	30181	6
14~16mm, オレンジ、配列 6x6	30183	2
17~20mm, 白、配列 4x5	30185	4
21~25mm、 青、配列 4x4	30187	2
26~30mm、 緑、配列 3x3	30189	4
1.5ml 微量遠心機用、 青、配列 4x6	30191	3
フルサイズ		
10~13mm、配列 6x12	30180BI	3
14~16mm、配列 6x12	30182	2
17~20mm、配列 4x10	30184	2
21~25mm、配列 4x10	30186	1
26~30mm、配列 3x8	30188	2
1.5ml 微量遠心機用、 配列 8x12	30190	2

<u>Solaris 2000 I / 2000 Rプラットフォーム用</u>

表 18: Solaris 2000 I / 2000 Rプラットフォーム用試験管立て製品

	製品番号	標準
		18 x 18
ハーフサイズ		
10~13mm, 赤、配列 6x6	30181	12
14~16mm, オレンジ、配列 6x6	30183	8
17~20mm, 白、配列 4x5	30185	10
21~25mm、 青、配列 4x4	30187	8
26~30mm、 緑、配列 3x3	30189	9
1.5ml 微量遠心機用、 青、配列 4x6	30191	8
フルサイズ		
10~13mm、配列 6x12	30180BI	7
14~16mm、配列 6x12	30182	4
17~20mm、配列 4x10	30184	5
21~25mm、配列 4x10	30186	3
26~30mm、配列 3x8	30188	4
1.5ml 微量遠心機用、 配列 8x12	30190	5

<u>Solaris 4000 I / 4000 Rプラットフォーム用</u>

表 19: Solaris 4000 I / 4000 Rプラットフォーム用試験管立て製品

1. 2. 5. 角度調節式試験管立てホルダー

<u>Solaris 2000プラットフォーム用</u>



	製品番号		標準		デュアル	スタック
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
10~13mm、72か所	236090	3	4	6	6	8
16~20mm、40か所	236091	3	4	6	5	8
21~25mm、40か所	236092	3	3	4	5	6
26~30mm、24か所	236093	3	3	4	5	6

表 20: Solaris 2000プラットフォーム用角度調節式試験管立てホルダー製品

<u>Solaris 4000プラットフォーム用</u>

	製品番号	標	뾑	デュアルスタック	
		18 x 30	36 x 24	18 x 30	
10~13mm、72か所	236090	7	14	14	
16~20mm、40か所	236091	7	14	14	
21~25mm、40か所	236092	5	9	10	
26~30mm、24か所	236093	6	9	12	

表 21: Solaris 4000プラットフォーム用角度調節式試験管立てホルダー製品

<u>Solaris 2000 I / 2000 Rプラットフォーム用</u>

	製品番号	標準
		11 x 14
10~13mm、72か所	236090	2
16~20mm、40か所	236091	2
21~25mm、40か所	236092	1
26~30mm、24か所	236093	1

表 22: Solaris 2000 I/2000 Rプラットフォーム用角度調節式試験管立てホルダー製品

<u>Solaris 4000 I / 4000 Rプラットフォーム用</u>

	製品番号	標準
		18 x 18
10~13mm、72か所	236090	4
16~20mm、40か所	236091	3
21~25mm、40か所	236092	3
26~30mm、24か所	236093	3

表 23: Solaris 4000 I/4000 Rプラットフォーム用角度調節式試験管立てホルダー製品

1. 2. 6. 調節式容器クランプ

<u>Solaris 2000プラットフォーム用</u>



	製品番号 標準			デュアル	スタック	
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
調節式一段式容器クランプ 11x14インチ(約30x36cm)	75004104	-	-	-	-	-
調節式二段式容器クランプ 12x14インチ(約30x36cm)	75004102	1	1	1	-	2
調節式二段式容器クランプ 18x18インチ(約46x46cm)	75004103	-	1	1	-	-
調節式一段式容器クランプ	75004101	4	4	6	6	8

表 24: Solaris 2000プラットフォーム用調節式容器クランプ製品

<u>Solaris 4000プラットフォーム用</u>

	製品番号	標準		デュアルスタック	
		18 x 30	36 x 24	18 x 30	
調節式一段式容器クランプ 11x14インチ(約30x36cm)	75004104	-	-	-	
調節式二段式容器クランプ 12x14インチ(約30x36cm)	75004102	2	2	4	
調節式二段式容器クランプ 18x18インチ(約46x46cm)	75004103	1	2	2	
調節式一段式容器クランプ	75004101	8	12	16	

表 25: Solaris 4000プラットフォーム用調節式容器クランプ製品

<u>Solaris 2000 I / 2000 Rプラットフォーム用</u>

	製品番号	標準
		11 x 14
調節式一段式容器クランプ 11x14インチ(約30x36cm)	75004104	1
調節式二段式容器クランプ 12x14インチ(約30x36cm)	75004102	-
調節式二段式容器クランプ 18x18インチ(約46x46cm)	75004103	-
調節式一段式容器クランプ	75004101	2

表 26: Solaris 2000 I/2000 Rプラットフォーム用調節式容器クランプ

<u>Solaris 4000 I / 4000 Rプラットフォーム用</u>

	製品番号	標準
		18 x 18
調節式一段式容器クランプ 11x14インチ(約30x36cm)	75004104	-
調節式二段式容器クランプ 12x14インチ(約30x36cm)	75004102	1
調節式二段式容器クランプ 18x18インチ(約46x46cm)	75004103	1
調節式一段式容器クランプ	75004101	4

表 27: Solaris 4000 I/4000 Rプラットフォーム用調節式容器クランプ

1. 2. 7. 角型メディアクランプ

<u>Solaris 2000プラットフォーム用</u>



	製品番号		標準		デュアル	スタック
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
125ml 角型ボトルクランプ	75004106	12	16	24	21	32
250ml 角型ボトルクランプ	75004107	9	16	22	18	32
500ml 角型ボトルクランプ	75004108	6	9	12	11	18
1000ml 角型ボトルクランプ	75004109	4	9	12	8	18

表 28: Solaris 2000プラットフォーム用角型メディアクランプ製品

<u>Solaris 4000プラットフォーム用</u>

	製品番号	標準		デュアルスタック	
		18 x 30	36 x 24	18 x 30	
125ml 角型ボトルクランプ	75004106	28	54	56	
250ml 角型ボトルクランプ	75004107	28	52	56	
500ml 角型ボトルクランプ	75004108	15	24	30	
1000ml 角型ボトルクランプ	75004109	15	24	28	

表 29: Solaris 4000プラットフォーム用角型メディアクランプ製品

<u>Solaris 2000 I / 2000 Rプラットフォーム用</u>

	製品番号	標準
		11 x 14
125ml 角型ボトルクランプ	75004106	9
250ml 角型ボトルクランプ	75004107	7
500ml 角型ボトルクランプ	75004108	3
1000ml 角型ボトルクランプ	75004109	2

表 30: Solaris 2000 I/2000 Rプラットフォーム用角型メディアクランプ製品

<u>Solaris 4000 I / 4000 Rプラットフォーム用</u>

	製品番号	標準
		18 x 18
125ml 角型ボトルクランプ	75004106	16
250ml 角型ボトルクランプ	75004107	16
500ml 角型ボトルクランプ	75004108	9
1000ml 角型ボトルクランプ	75004109	9

表 31: Solaris 4000 I/4000 Rプラットフォーム用角型メディアクランプ製品

1. 2. 8. Nalgeneビーカーラック

<u>Solaris 2000プラットフォーム用</u>



	製品番号		標準		デュアル	スタック
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
寸法 11x14インチ (約30x36cm)/12x14インチ (約30x36cm)			Ę	ラック数/立本	数	
Nalgene 30mlビーカー、 立本数36本	75004116	1 / 36	-	-	2 / 72	-
Nalgene 50mlビーカー、 立本数28本	75004129	1 / 28	-	-	2 / 56	-
Nalgene 100mlビーカー、 立本数16本	75004118	1 / 16	-	-	2 / 32	-
Nalgene 250mlビーカー、 立本数9本	75004119	1 / 9	-	-	2 / 18	-
Nalgene 400mlビーカー、 立本数8本	75004120	1 / 8	-	-	2 / 16	-
Nalgene 600mlビーカー、 立本数5本	75004121	1 / 5	-	-	2 / 10	-
寸法 9x18インチ (約23x46cm)			Ę	ラック数/立本	数	
Nalgene 30mlビーカー、 立本数32本	75004110	-	2 / 64	2 / 64	-	4 / 128
Nalgene 50mlビーカー、 立本数24本	75004128	-	2 / 48	2 / 48	-	4 / 96
Nalgene 100mlビーカー、 立本数15本	75004112	-	2 / 30	2 / 30	-	4 / 60
Nalgene 250mlビーカー、 立本数8本	75004113	-	2 / 16	2 / 16	-	4 / 32
Nalgene 400mlビーカー、 立本数6本	75004114	-	2 / 12	2 / 12	-	4 / 24
Nalgene 600mlビーカー、 立本数4本	75004115	-	2 / 8	2 / 8	-	4 / 16

表 32: Solaris 2000プラットフォーム用Nalgeneビーカーラック製品

34 / 138

<u>Solaris 4000プラットフォーム用</u>

	製品番号	標準		デュアルスタック
		18 x 30	36 x 24	18 x 30
寸法 11x14インチ (約30x36cm)/12x14インチ (約30x36cm)			ラック数/立本	≍数
Nalgene 30mlビーカー、 立本数36本	75004116	-	-	-
Nalgene 50mlビーカー、 立本数28本	75004129	-	-	-
Nalgene 100mlビーカー、 立本数15本	75004118	-	-	-
Nalgene 250mlビーカー、 立本数9本	75004119	-	-	-
Nalgene 400mlビーカー、 立本数8本	75004120	-	-	-
Nalgene 600mlビーカー、 立本数5本	75004121	-	-	-
寸法 9x18インチ(約 23x46cm)			ラック数/立本	蒸数
Nalgene 30mlビーカー、 立本数32本	75004110	3 / 96	4 / 128	6 / 192
Nalgene 50mlビーカー、 立本数24本	75004128	3 / 72	4 / 96	6 / 144
Nalgene 100mlビーカー、 立本数15本	75004112	3 / 45	4 / 60	6 / 90
Nalgene 250mlビーカー、 立本数8本	75004113	3 / 24	4 / 32	6 / 48
Nalgene 400mlビーカー、 立本数6本	75004114	3 / 18	4 / 24	6 / 36
Nalgene 600mlビーカー、 立本数4本	75004115	3 / 12	4 / 16	6 / 24

表 33: Solaris 4000プラットフォーム用Nalgeneビーカーラック製品

35 / 138

<u>Solaris 2000 I / 2000 Rプラットフォーム用</u>

	製品番号	標準
		11 x 14
寸法 11x14インチ(約30x36cm)/12x14 インチ(約30x36cm)		ラック数/立本数
Nalgene 30mlビーカー、立本数36本	75004116	1 / 36
Nalgene 50mlビーカー、立本数28本	75004129	1 / 28
Nalgene 100mlビーカー、立本数16本	75004118	1 / 16
Nalgene 250mlビーカー、立本数9本	75004119	1 / 9
Nalgene 400mlビーカー、立本数8本	75004120	1 / 8
Nalgene 600mlビーカー、立本数5本	75004121	1 / 5
寸法 9x18インチ (約23x46cm)		ラック数/立本数
Nalgene 30mlビーカー、立本数32本	75004110	-
Nalgene 50mlビーカー、立本数24本	75004128	-
Nalgene 100mlビーカー、立本数15本	75004112	-
Nalgene 250mlビーカー、立本数8本	75004113	-
Nalgene 400mlビーカー、立本数6本	75004114	-
Nalgene 600mlビーカー、立本数4本	75004115	-

表 34: Solaris 2000 I/2000 Rプラットフォーム用Nalgeneビーカーラック製品
付属品

<u>Solaris 4000 I / 4000 Rプラットフォーム用</u>

	製品番号	標準
		18 x 18
寸法 11x14インチ(約30x36cm)/12x14 インチ(約30x36cm)		ラック数/立本数
Nalgene 30mlビーカー、立本数36本	75004116	-
Nalgene 50mlビーカー、立本数28本	75004129	-
Nalgene 100mlビーカー、立本数16本	75004118	-
Nalgene 250mlビーカー、立本数9本	75004119	-
Nalgene 400mlビーカー、立本数8本	75004120	-
Nalgene 600mlビーカー、立本数5本	75004121	-
寸法 9x18インチ (約23x46cm)		ラック数/立本数
Nalgene 30mlビーカー、立本数32本	75004110	2 / 64
Nalgene 50mlビーカー、立本数24本	75004128	2 / 48
Nalgene 100mlビーカー、立本数15本	75004112	2 / 30
Nalgene 250mlビーカー、立本数8本	75004113	2 / 16
Nalgene 400mlビーカー、立本数6本	75004114	2 / 12
Nalgene 600mlビーカー、立本数4本	75004115	2 / 8

表 35: Solaris 4000 I/4000 Rプラットフォーム用Nalgeneビーカーラック製品

1. 2. 9. 分液ロートクランプ

<u>Solaris 2000プラットフォーム用</u>



	製品番号		標準		デュアル	スタック
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
分液ロート 250ml~2,000ml 用シングルクランプ	75004125	2	3	4	2	5
分液ロート垂直取付用ク ランプ		4	4	6	-	-

表 36: Solaris 2000プラットフォーム用分液ロートクランプ製品

<u>Solaris 4000プラットフォーム用</u>

	製品番号	標準		デュアルスタック
		18 x 30	36 x 24	18 x 30
分液ロート 250ml~2,000ml 用シングルクランプ	75004125	5	7	9
分液ロート垂直取付用クランプ		8	12	-

表 37: Solaris 4000プラットフォーム用分液ロートクランプ製品

<u>Solaris 2000 I / 2000 Rプラットフォーム用</u>

	製品番号	標準
		11 x 14
分液ロート 250ml~2,000ml 用シングルクランプ	75004125	
分液ロート垂直取付用クランプ		-

表 38: Solaris 2000 I/2000 Rプラットフォーム用分液ロートクランプ製品

<u>Solaris 4000 I / 4000 Rプラットフォーム用</u>

	製品番号	標準
		18 x 18
分液ロート 250ml~2,000ml 用シングルクランプ	75004125	-
分液ロート垂直取付用クランプ		-

表 39: Solaris 4000 I/4000 Rプラットフォーム用分液ロートクランプ製品

38 / 138

1. 2. 10. ユーティリティトレイ

<u>Solaris 2000プラットフォーム用</u>

www.

	製品番号	標準		デュアルスタック		
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
11x14インチ (約30x36cm)/12x14 インチ(約30x36cm) ユーティリティトレイ	75004123	1	-	-	2	-
9x18インチ(約23x46cm) ユーティリティトレイ	75004122	-	2	3	-	4

表 40: Solaris 2000プラットフォーム用ユーティリティトレイ製品

<u>Solaris 4000プラットフォーム用</u>

	製品番号	標準		デュアルスタック
		18 x 30	36 x 24	18 x 30
11x14インチ (約30x36cm)/12x14インチ (約30x36cm)ユーティリテ ィトレイ	75004123	-	-	-
9x18インチ(約23x46cm)ユーテ ィリティトレイ	75004122	3	6	6

表 41: Solaris 4000プラットフォーム用ユーティリティトレイ製品

<u>Solaris 2000 I / 2000 Rプラットフォーム用</u>

	製品番号	標準
		11 x 14
11x14インチ(約30x36cm)/12x14インチ (約30x36cm)ユーティリティトレイ	75004123	1
9x18インチ(約23x46cm)ユーティリティトレイ	75004122	-

表 42: Solaris 2000 I/2000 Rプラットフォーム用ユーティリティトレイ製品

<u>Solaris 4000 I / 4000 Rプラットフォーム用</u>

	製品番号	標準
		18 x 18
11x14インチ(約30x36cm)/12x14インチ (約30x36cm)ユーティリティトレイ	75004123	-
9x18インチ (約23x46cm) ユーティリティトレイ	75004122	2

表 43: Solaris 4000 I/4000 Rプラットフォーム用ユーティリティトレイ製品

39 / 138

ユーティリティトレイ付属品

	製品番号	標準
0リングのスペアセット (ユーティリティトレイ)	75004132	-

 \checkmark

表 44: ユーティリティトレイ付属品

1. 2. 11. 粘着マット

粘着マットの選択に関する一般的な注意事項

- 寸法170x280mmのマットは、11x14インチおよび12x14インチのプ ラットフォーム用にご利用ください。
- 寸法200x200mmのマットは、より大きな18x18インチ、18x24インチ、 30x18インチ、36x24インチのプラットフォーム用にご利用ください。
- マットは適切なサイズに切って使用することができます。

Solaris 2000プラットフォーム用

	製品番号		標準		デュアル ク樽	レスタッ 標準
		12 x 14	18 x 18	18 x 24	12 x 14	18 x 18
200x200mm 高粘着マット、2枚 入パック	75004126	1	4	6	2	8
280x170mm 高粘着マット、2枚 入パック	75004127	2	2	4	4	4
200x200mm 低粘着マット、2枚 入パック	75004111	1	4	6	2	8
280x170mm 低粘着マット、2枚 入パック	75004117	2	2	4	4	4

マットは2枚入パックで販売されています。このパックは2枚入ですが、プラットフォームに敷くマットの枚数は1枚です。

表 45: Solaris 2000プラットフォーム用粘着マット製品

<u>Solaris 4000プラットフォーム用</u>

		製品番号	標準		デュアルスタ ック
			8 x 30	36 x 24	18 x 30
	200x200mm 高粘着マット、2枚入パック	75004126	6	12	12
	280x170mm 高粘着マット、2枚入パック	75004127	6	9	12
	200x200mm 低粘着マット、2枚入パック	75004111	6	12	12
	280x170mm 低粘着マット、2枚入パック	75004117	6	9	12
_				*	

マットは2枚入パックで販売されています。このパックは2枚入ですが、プラットフォームに敷くマットの枚数は1枚です。

表 46: Solaris 4000プラットフォーム用粘着マット製品

<u>Solaris 2000 I / 2000 Rプラットフォーム用</u>

	製品番号	標準
		11 x 14
200x200mm 高粘着マット、2枚入パック	75004126	1
280x170mm 高粘着マット、2枚入パック	75004127	2
200x200mm 低粘着マット、2枚入パック	75004111	1
280x170mm 低粘着マット、2枚入パック	75004117	2

マットは2枚入バックで販売されています。このバックは2枚入ですが、ブラットフォームに敷くマットの枚数は1枚です。 表 47: Solaris 2000 I/2000 Rプラットフォーム用粘着マット製品

<u>Solaris 4000 I / 4000 Rプラットフォーム用</u>

	製品番号	標準
		18 x 18
200x200mm 高粘着マット、2枚入パック	75004126	4
280x170mm 高粘着マット、2枚入パック	75004127	2
200x200mm 低粘着マット、2枚入パック	75004111	4
280x170mm 低粘着マット、2枚入パック	75004117	2

マットは2枚入パックで販売されています。このパックは2枚入ですが、プラットフォームに敷くマットの枚数は1枚です。

表 48: Solaris 4000 I/4000 Rプラットフォーム用粘着マット製品

41 / 138

1. 2. 12. その他の付属品

1. 2. 12. その他の付属		and the second s
	製品番号	
粘着マット用ローリングアプリケーター	75004124	
付属品用ドライバー	75004131	
ガス供給マニホールド (Solaris 2000 I/2000 R用)	SK2000-8GM	
ガス供給マニホールド (Solaris 4000 I/4000 R用)	SK4000-8GM	
クランプスペアキット(ネジ)	SK0010	
Solaris 2000、2000 I/R、4000 I/R用のプ ラットフォームスペアキット(プラットフ ォーム用ネジ、工具)	SK0100	
SK4000用プラットフォームスペアキット (プラットフォーム用ネジ、工具)	SK0101	

表 49: その他の付属品

1. 3. 指令および 規格

地域	指令	規格
ヨーロッパ	2006/42/EC	EN 61010-1
	Machinery Directive	IEC 61010-2-051
		EN 61326-1 クラスB
	2011/65/EU RoHS	EN ISO 14971
	電気・電子機器に含まれる特定	EN ISO 9001
	有吉物貨の使用制限 - 関9 る指 今	<u>Solaris 2000 I/2000 R/</u>
		<u>4000 I/4000 Rのみ対象</u>
	<u>休啶口标·</u> 2014/25/EU 低電口	IEC 61010-2-010
	2014/33/L0 低电圧	
	2014/30/20 电燃间立住 (2007	ANSI/III 61010-1
山木		ANSI/ 02 01010 1
		CFR 4/ FCC 15 EMC
		EN ISO 14971
		EN ISO 9001
		<u>Solaris 2000 I/2000 R/</u>
		<u>4000 I/4000 Rのみ対象</u>
口平		
		1EC 61010-2-051
		IEC 61326-1 クラスB
		EN ISO 14971
		EN ISO 9001
		<u>Solaris 2000 I/2000 R/</u>
		<u>4000 I/4000 Rのみ対象</u> IEC 61010-2-010

表 50: 指令および規格

留意: 本機器は、FCC規定の第15章に定められたクラスAデジタル機器に関する規制要件に 基づいて所定の試験が実施され、これに適合するものと認定されています。この規制は、 商用環境下での設置に際し、有害な電波障害に対する合理的な保護を規定するために策定 されました。本機器は高周波エネルギーを発生させ、利用するもので、かつ外部に放射す ることがあるので、取扱説明書に従って設置、使用されない場合は、無線通信に有害な支 障を引き起こす可能性があります。そのため、住宅領域でこの機器を使用すると有害な電 波障害が生じる可能性があり、その場合にはユーザーが自己負担にてその障害の修復をそ の障害の修復をすることになります。

1. 4. 製品概要

1. 4. 1. Solaris 2000



プラットフォーム取付板;② グラフ表示のあるユーザーインターフェース;
 電源スイッチ;④ 電源;⑤ ヒューズ;⑥ イーサネットポート;⑦ USBポート
 Solaris 2000概要

1. 4. 2. Solaris 4000



 プラットフォーム取付板:② グラフ表示のあるユーザーインターフェース:③ 電源 スイッチ:④ 電源:⑤ ヒューズ(リセット機能付):⑥ イーサネットポート:⑦ USB ポート

図 2: Solaris 4000概要

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



① プラットフォーム取付板:② グラフ表示のあるユーザーインターフェース:③ 電源スイ ッチ:④ アクセスポート:⑤ 電源接続部:⑥ ヒューズ(リセット機能付):⑦ イーサネ ットポート:⑧ USBポート:⑨ フード

図 3: Solaris 2000 I/2000 R概観

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



① プラットフォーム取付板; ② グラフ表示のあるユーザーインターフェース; ③ 電源ス イッチ; ④ アクセスポート; ⑤ 電源接続部; ⑥ ヒューズ(リセット機能付); ⑦ イーサ ネットポート; ⑧ USBポート; ⑨ フード

図 4: Solaris 4000 I/4000 R概観

1. 4. 5. 接続

開放型シェーカー



① 電源: ② ヒューズ: ③ イーサネットポート

図 5: 背面図 (左: Solaris 4000、右: Solaris 2000)

恒温シェーカー(加熱式)および冷却シェーカー(冷却・加熱式)



① 電源:
 ② ヒューズ:
 ③ イーサネットポート:
 ④ ドレンチューブ
 図 6: 背面図(左: Solaris 2000 I / 2000 R, 右: Solaris 4000 I / 4000 R)

電源

シェーカーには、各シェーカーの仕様に対応する電源が必要です。電源ケーブル は同梱されています。

製品番号	シェーカー	仕様
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

図 7: シェーカー別電源仕様

警告 不適切な電源または電源プラグによる損傷。シェーカーを接続するコンセントが有効な電気機器の安全規格に従って適切に接地されていることを必ず確認してください。損傷した電源ケーブルや不適切な定格の電源ケーブルでシェーカーを接続しないでください。

コンセントプラグは、如何なる時もすぐに届くようにして下さい。

感電の危険性を排除するために、シェーカー周囲の表面が乾いていることを確認 してください。液体の流出・飛散事故が発生した場合は、シェーカーのプラグを 外し、流出した液体を清掃し、健康被害や生物学的危険性を排除してから続行し てください。

シェーカー不使用時はプラグを外してください。

ヒューズ

ヒューズは、過電流状態が発生したために切れたとき、戻すことができます。シ ェーカーはヒューズを定位置に戻すまで動作しません。

イーサネット

シェーカーには、ローカルエリアネットワーク(LAN)への接続に使用できる RJ45イーサネットポートが内蔵されています。接続する機器は、RJ45イーサネ ットポートが装備された規格IEC 60950-1に準拠したもののみを使用してくださ い。RJ45イーサネットポートは、今後、ソフトウェアの適切なアップデートが利 用可能になったとき使用するために装備されています。

アクセスポート

温度調節機能付シェーカーのモデルには、アクセスポートと呼ばれる穴が側面に 2つあります。アクセスポートは通常、キャップで塞がれています。キャップを 取り外して、予備の温度センサーなどの追加の器具を取り付けたり、ケーブルや ガス供給マニホールドのチューブなどのチューブ類を通したりできます。

47 / 138

USB

シェーカーにはUSB-A 2.0ポートが2個内蔵されており、市販のUSBドライブが使 用できます。接続する機器は、USBポートが装備された規格IEC 60950-1に準拠し たもののみを使用してください。

内部過熱防止策

温度調節機能付シェーカーのモデルには、発熱体の横に手動リセットのサーモス タットが取り付けられています。この温度監視装置により、空気循環用ファンが 故障した場合に、シェーカーのキャビネット内で温度が過昇しないようにヒータ ーが遮断されます。

手動リセットのサーモスタットでは、シェーカーを元の動作状態に戻すには、手 動でセットし直す必要があります。この手順は、Thermo Fisher Scientificカス タマーサービスだけが実行できます。 開梱

2. 運搬と設置

i

ユーザーは責任を持って、シェーカーが正しく設置されてい るか確認してください。

注

必ず納入時に輸送用カートンを検査してください。受け取ったら、開梱する前に 輸送中の損傷を慎重に点検してください。損傷が見つかった場合は、運送業者が 納品受領証の顧客用控えに損傷を明記し、署名することになっています。

カートンを慎重に開き、梱包資材を捨てる前にすべての部品(「表 51: 納品 内容」)が同梱されているか確認してください。開梱後、損傷が見つかった場合 は、運送業者に報告して損傷検査を依頼してください。

重要: 輸送品の受領後数日以内に損傷検査を要請しなかった場合は、損害に 対して運送業者を免責することになります。損傷検査は必ず要請してください。

2.1. 開梱

開梱時に梱包リストに記載してあるユニットー式を受け取ったか確認してください。すべて納得できるまで梱包材を捨てないでください。

納品内容

品目	数
シェーカー	1
電源コード	1
標準プラットフォーム	1*
プラットフォーム用ネジ	
- Solaris 4000	4*
- その他のモデル	3*
取扱説明書(印刷物、英語)	1
USBのマニュアル	1
ロックツール(「ハンドルレンチ)	1
* 2000 I、2000 R、4000 I、4000 Rシェーカーには	、シェーカーにプ
ラットフォームがあらかじめ組み込まれています。	プラットフォーム
やネジなどの個々の部品はありません。	

表 51: 納品内容

納品内容の物品が同梱されていない場合は、Thermo Fisher Scientificまでご連 絡ください。

2. 2. 設置場所

注意 プラスチックは紫外線にさらされると安定性が低下するため、損傷に対する防護が 衰える可能性があります。シェーカーやプラスチック製の付属品は直射日光や紫外線の光 源などに露出させないでください。

シェーカーは動作中に付属品やサンプルを含めたシェーカーの重量を支えること ができる水平な実験台やベンチなどに置いてください。シェーカーは銘板に記載 してある要件に合ったコンセントの近くに置いてください。装置を設置するとき は周囲の間隔を空けて、空気の自然対流を妨げることなく、付属品の取り付けが 可能でユーザーが使いやすいようにしてください。

装置を設置するには、以下の要件に注意してください。

- 開放型シェーカーを配置するときは、可動する部分の動きを妨げないように 十分な間隔を確保し、圧搾による負傷や隣接する機器の破損を防ぐために、 プラットフォームのすべての4辺から8cm(3インチ)間隔を空けます。開放 型シェーカーのプラットフォームが台の縁と重なる場合があります。
- シェーカーは振動を引き起こします。危険防止ゾーンには、敏感な機器、危険品または有害物質を保管しないでください。



図 8: 周囲8cm (3インチ)の間隔を空けた開放型シェーカーの配置方法

温度調節機能付シェーカーにおいて、適切な換気で性能を最大に発揮するには、 以下の間隔要件に従って設置してください。

- Solaris 2000 I・4000 Iなどの恒温シェーカー(加熱式)には、図 9の左図 に示すように、台の縁からすべての4辺に対して8cm(3インチ)の空間が必 要です。
- Solaris 2000 R・4000 Rなどの冷却シェーカー(冷却・加熱式)では、図 9 の左図に示すように、フードの上に何もない空間がある場合は、台の縁から すべての4辺に対して8cm(3インチ)間隔を空けて操作します。

 Solaris 2000 R・4000 Rなどの冷却シェーカー(冷却・加熱式)の上の空間 が遮られている場合は、図 9の右図に示すように、台の縁の周囲からすべて の4辺に対して30cm(12インチ)の空間に増やす必要があります。



図 9: 恒温シェーカー(加熱式)(左)と冷却シェーカー(冷却・加熱式)(右)の間隔を空けた設置図。冷却シェーカー(冷却・加熱式)の上に何もない場合は、左の画像が適用されます。

警告 振とう中は、有害物質を区域外に隔離してください。 支持面に必要なことは以下の通りです。

- » 清潔であること
- » 安定性が良く頑丈であり、固定されていて共振がないこと
- » シェーカーを水平に設置するのに最適であること
- » シェーカーの重量に耐えられること
- シェーカーは室内専用です。
- 高温になる場所や直射日光があたる場所を避けてください。
- 設置場所は、常時良く換気されている場所であること。
- コンセントプラグは、如何なる時もすぐに届くようにして下さい。

2.3. 運搬

注意 シェーカーの落下による身体的危害。シェーカーは必ず両側から持ち上げてください。シェーカーはフロントパネルや取り付けてあるプラットフォームを掴んで持ち上げないでください。温度調節機能付シェーカーは、透明なフードに付いているハンドルで持ち 上げないでください。

注意 台の縁の中で水が結露すると、シェーカーが安全に使用できません。これは、シェ ーカーが涼しい場所から湿度の高い暖かい場所に移動したときに発生します。たとえば、 保管場所や運搬トラック から実験室に移動した場合です。そのため、操作する前に少なく とも2時間乾燥させてください。

注 開放型シェーカーモデルでは、シェーカーを移動する前に、必ずプラットフォーム、 その上の積載物、すべての付属品を取り外してください。プラットフォームを取り外さな いと、プラットフォーム取付板や振動機構が損傷する可能性があります。温度調節機能付 シェーカーでは、プラットフォームから積載物を取り除くだけで十分です。

注 ユニットへの衝撃による破損の危険性あり。シェーカーは直立させた状態で、できれ ば専用の梱包にて輸送してください。

シェーカーの取扱い

シェーカーを取り扱うときは、以下の事柄を確認してください。

- シェーカーは左右両側面で持ち上げます。正面と背面で持ち上げないこと。
- シェーカーの重量(11ページの「技術仕様」参照)に応じて、できるだけ 多くの人で運ぶように割り当てます。1人で運ばないこと。







図 10: シェーカーの持ち運び方

警告 シェーカーは必ず両側から持ち上げてください。シェーカーを正面や背面から持ち 上げないでください。シェーカーはかなり重量があります(11ページの「技術仕様」参 照)。シェーカーを持ち上げて運ぶときは、最低2人を割り当ててください。

注意 温度調節機能付シェーカーは、透明なフードに付いているハンドルをつかんで持ち 上げないでください。ハンドルが破損し、シェーカーが地面に落ちて、重傷を負う危険性 があります。

2. 4. 水平調整方法

シェーカーには本体を水平にする機能がありません。シェーカーを適切に設置す るために水平な支持面が必要です。

注 シェーカーを水平に調整するために、シェーカーの脚と支持面の間にくさびなどの平 らなものを置かないでください。

2.5. 電源

シェーカーには、各シェーカーの仕様に対応する電源が必要です。電源ケーブル は同梱されています。詳細については、47ページの図 7ページの図7を確認し てください。

警告 不適切な電源または電源プラグによる損傷。シェーカーが適切に接地されたソケットに差し込まれていることを確認してください。損傷した電源ケーブルや不適切な定格の電源ケーブルでシェーカーを接続しないでください。

注 電磁放射は、ディスプレイに電波干渉を引き起こす可能性があります。ただし、これ によってデバイスの破損、機能の制限や変更が引き起こされることはありません。電磁放 射による電波干渉を避けるために、携帯電話などのモバイルデバイスを本機の近辺に持ち 込まないでください。他の高電力の機器と一緒の回路上で本機を動作しないでください。 同じ電源タップで複数の機器を動作しないでください。

シェーカーを電源に接続するには、次の手順で行います。

- 1. 本体右側にある電源スイッチを切ります。
- ケーブルの仕様が、ご利用の国での安全基準に沿っているかを確認してくだ さい。
- 3. 電圧と周波数が、定格板に記載のある数字と同様かを確認してください。

コンセントプラグは、如何なる時もすぐに届くようにして下さい。

感電の危険性を排除するために、シェーカー周囲の表面が乾いていることを確認 してください。液体の流出・飛散事故が発生した場合は、シェーカーのプラグを 外し、流出した液体を清掃してから続行してください。

シェーカー不使用時はプラグを外してください。

53 / 138

2. 6. 初回起動

装置を使用する前に、以下の手順を行ってください。

 装置の電源を入れると、thermo scienceのロゴが表示されます。設定を開始 するをタップします。

ther	moscientific	
	Start Setup	

図 11: 初回起動プロンプト

2. 言語画面で使用する言語を選択します。次へをタップします。

Language			
Deutsch			
English			
Fspañol			
Back Next			

図 12: 初回起動 - 言語

 オプションとして、ユニット名をダイアログボックスの「ユニット名」の欄 に入力します。次へをタップします。

		Unit M	Name	
	Unit Name:	Tap to enter		
		Back	Next	
义	13: 初回	起動	ユニット名	

 オプションとして、テキストフィールドに都市と国を入力します。3文字を 入力するとすぐに、一致する候補のリストが「都市と国」テキストフィール ドに表示されます。必要に応じて、1つを選択し、次へをタップします。

Set Region		
	Enter the city and country closest to your time zone	
City, Country: Tap to enter		
В	ack Next	

図 14: 初回起動 - 地域

5. 使用する日付フォーマットを選択します。次へをタップします。

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

図 15: 初回起動 - 日付表示フォーマットの設定

 ホイールピッカーの各ダイヤルを回して、現在の日付を選択します。次へを タップします。



図 16: 初回起動 - 日付の設定

 使用する時刻フォーマットを選択し、ホイールピッカーの各ダイヤルを回し て現在の時刻を設定します。次へをタップします。



図 17: 初回起動 - 時間の設定

恒温シェーカーおよび冷却シェーカー用: 使用する温度単位を選択します。次へをタップします。

Temperate	ure Units
	°C
0	°F
0	
Back	Next

図 18: 初回起動 - 温度単位の選択

 恒温シェーカーおよび冷却シェーカー用: 青色の矢印をクリックして、高 温アラームと緊急運転停止の閾値を設定します。次へをタップします。

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

- 図 19: 初回起動 高温・低温アラームの閾値の設定
- 10. 上記の手順で、低温アラームの閾値を設定します。次へをタップして続行します。

 管理者パスコードを設定せずに続行する場合は、オープンモードを選択した まま、次へをタップして続行します。

ここで管理者パスコードを設定する場合は、セキュアモードをタップします。

		,	Acce	ss Mode				
	۲	Open Mode Select to allow full a	ccess	to all functionality for all users.				
	Secure Mode Select to require users to enter the admin passcode to make changes to stored parameters, programs, and settings. (All users still able to use programs and set new run parameters.)							
		Back		Next				
义	20:	初回起動	_	アクセスモードの選択				

12. パスコードプロンプトが表示されたら、キーパッドを使用して管理者パスコ ードを入力し、次へをタップします。

	Admin F	Passcode
The unit h like to chan	as shipped with a s ge the standard adn passco	tandard passcode. If you would nin passcode, please enter a new de below.
If you do no fi	t wish to change the elds below blank ar	e passcode at this time, leave the d tap the Next button.

図 21: 初回起動 - アクセスモードの選択

- 次のパスコードプロンプトが表示されたら、管理者パスコードを入力します。
- 3番目のパスコードプロンプトが表示されたら、管理パスコードをもう一度 入力して確認します。

	Admin Passcode					Adm	in Passo	ode	
If you do not w field	ish to change Is below blan	e the passe k and tap t	code at this he Next bu	s time, leave the itton.	P	lease re-ent	er the ne	w passcode.	
*	*	*	*	*	*	*	*	*	*
	7	8	9			7	8	9	
	4	5	6			4	5	6	
	1	2	3			1	2	3	
	0	Clear				0	Clea	ar	

図 22: 初回起動 - 新規管理者パスコードの入力と確認

57 / 138

15. 確認画面が表示されたら、次へをタップして続行します。



- 16. シェーカーの一般的な物理的設置ガイドに関する画面が表示されます。次へ をタップします。

		Installation	Instructions
	SI s refe	hakers must be instal surface with sufficien rence the product ma proper installa	ied on a stable and level t air circulation. Please nual for instruction on the ion of this unit.
		Back	Next
义	24:	初回起動 -	設置ガイド

17. 「設定完了」ウィンドウが表示されます。次へをタップして終了します。

Setup C	omplete
Setup is o	complete.
We recommend calibra temperature for your app calibrate the ter	ting the unit at the ideal lication. Would you like to nperature now?
Calibration can always be accessed	ed under the main settings screen.
Calibrate	Finish

図 25: 初回起動 - 校正(較正)の設定または終了の設定

注 恒温シェーカー(加熱式)および冷却シェーカー(冷却・加熱式)では、校正を タップすると、126ページの「4.6.温度校正」の章で説明されている手順が実行で きます。

2. 7. 保管



シェーカーと付属品をしばらく使用しないときは、装置全体 を清掃し、必要に応じて消毒・除染を行ってください。シェ ーカーや付属品を不明確な汚染状態で放置しないでくださ

注意 い。手順に関してご不明な点がありましたら、Thermo Fisher Scientificカスタマーサービスまでお問い合わせください(123 ページの「クリーニング」、124ページの「消毒」、124ペー ジの「除染」)。

- シェーカーと付属品を保管する場合は清掃し、必要に応じて消毒・除染を行ってから保管してください。
- シェーカーと付属品は完全に乾かしてから保管してください。
- シェーカーは清潔でほこりのない場所に保管してください。
- シェーカーは倒れないように置いてください。
- シェーカーは直射日光の当たる場所に保管しないでください。

2. 8. 運送



注意

シェーカーと付属品を発送する場合は、装置全体のクリーニン グを行い、必要に応じて、殺菌または除染してから発送してく ださい。シェーカーや付属品を不明確な汚染状態で放置しない でください。手順に関してご不明な点がありましたら、Thermo Fisher Scientificカスタマーサービスまでお問い合わせくだ さい(123ページの「クリーニング」、124ページの「消毒」、

124ページの「除染」)。

シェーカーを発送するときは、以下の事柄を確認してください。

- シェーカーが清潔で除染されている必要があります。
- 除染証明で除染したことを裏付けること。除染証明書はThermo Fisher Scientificカスタマーサービスで取得できます。

3. 運転

3. 1. 電源のオン/オフ

本体右側にある電源スイッチを押して、シェーカーの電源をオン(I)またはオ フ(0)にします。

起動時に、タッチスクリーンにThermo Scientificのロゴが表示されます。

使用準備が完了するとタッチスクリーンにシェーカーの現在のステータスが表示 されます。

3. 2. グラフ表示のあるユーザーインターフェース

ホーム画面は、シェーカーのグラフィカルユーザーインターフェイス(GUI)に おけるデフォルト画面です。この画面はシェーカーを操作するための出発点にな ります。

ホーム画面では、以下の操作ができます。

- 振とう速度、振とう時間、振とう温度などの基本的な操作パラメーターの設定(温度調節機能付シェーカーのみ)
- シェーカーの開始および停止
- ステータス情報の表示およびアラームとアラートの管理

 ステータス情報と設定オプションを含む他の画面へのナビゲート 画面の内容は、開放型バージョンと温度調節付バージョンでわずかに異なります。

開放型シェーカーのホーム画面は、図 26のようになります。



図 26: 開放型シェーカーのGUIホーム画面

温度調節機能付シェーカーのホーム画面には、図 27のように、温度調節に関す る表示が追加されています。



図 27: 温度調節機能付シェーカーのGUIホーム画面

3. 2. 1. シェーカーの基本操作パラメーターの設定

ホーム画面で1回タップするだけで、主要なシェーカーの操作パラメーターの設 定画面が表示されます。

設定速度

実際速度フィールド(図 26の⑥または図 27の⑦)の任意の場所をタップして、図 28に示す設定速度画面を開きます。



図 28: 設定速度

- ホイールピッカーの各ダイヤルの上または下の矢印をタップして、目的とする速度を設定します。または、ホイールピッカーの各ダイヤルを個別に回転して、目的とする設定速度を変更することもできます。
- 3. 保存をタップします。

注 入力した設定速度が範囲外であることを通知するポップアップ画面が表示された 場合は、11ページの「仕様」の章に明記されているシェーカーの対応動作速度の範 囲外になります。設定を修正して続行します。

 メイン画面に戻るために表示される「設定が保存されました」確認ダイアロ グを了解します。

軌道計算機:従来の機器設定からおおよその速度の算出

「速度設定値」ダイアログ内から、速度を計算をタップして軌道計算機を開くこ ともできます。軌道計算機は、異なる機械設計のシェーカーから、本書で説明さ れているSolarisシリーズのシェーカーに移行するときに役立ちます。この計算 機では、軌道サイズの違いを用いて、同様の結果を得るのに必要な設定速度を概 算することができます。

注 軌道計算機で得られた結果を「そののまま」生産試料に使用するのではなく、一連の 試験を実施して確認してください。細胞は、異なる軌道に移動したときに物理現象の変化 につれて、せん断応力に応じて、より速くまたはより遅く成長したり、異なるタンパク質 を発現したり、損傷(または死滅)したりする可能性があります。

1. 測定単位をミリメートルまたはインチで選択します。



図 29: 軌道計算機

グラフ表示のあるユーザーインターフェース

 前回の軌道サイズを選択します。計算をタップします。速度の計算中は次の 画面が表示されます。



図 30: 軌道計算の結果

- 3. OKをタップします。
- また、推奨速度は軌道サイズをカスタマイズしても計算できます。カスタム 軌道を計算するには、「その他」を選択します。
- 5. 軌道サイズを入力します。



図 31: カスタマイズした軌道サイズの計算

6. 計算をタップします。

「速度計算」ウィンドウに、入力したサイズの速度が表示されます。

ランタイム(運転時間)の設定

シェーカーは連続モードまたはタイマーモードのどちらかで実行できます。連続 モードの場合は、ユーザーの判断でシェーカーを手動で停止します。タイマーモ ードの場合は、タイマーが切れるときにシェーカーの駆動が自動的に停止しま す。タイマーの表示について、以下のような選択できます。

- 経過時間: スタートボタンを押してからシェーカーが稼働している時間
- 残り時間: タイマーが切れるまでシェーカーが動作し続ける時間
- 時間フィールド(図 26および図 27の⑤)をタップして、図 32に示す時間モ ード画面を開きます。

2. タイマーを選択します。



図 32: 時間モード

- 時間の設定ボックスにある時間と分のに数字を入力します。
- 4. 保存をタップします。
- メイン画面に戻るには、表示される「設定が保存されました」確認ダイアロ グを了解します。

温度設定

温度調節付モデルの場合は、温度設定値フィールドで、いつでもアプリケーショ ンの温度を設定できます。

- 温度設定値 フィールド(図 27の項目⑥の一番下の欄)の任意の場所をタッ プして、図 33に示す温度設定値画面を開きます。
- シェーカーが維持する温度を決めて、ホイールピッカーの各ダイヤルの 上または下の矢印をタップし、デフォルト温度を事前に設定します。 または、ホイールピッカーの各ダイヤルを個別に回転して、目的とする温度 を変更することもできます。



図 33: 温度設定値の設定: 読み出し校正の表示がある画面

3. 保存をタップします。

注 入力した設定温度が範囲外であることを通知するポップアップ画面が表示された 場合は、11ページの「仕様」の章に明記されているシェーカーの対応動作温度の範 囲外になります。設定を修正して続行します。

注 ご利用時の周囲温度が原因で、選択した温度が範囲外であることを通知するポッ プアップ画面が表示された場合は、11ページの「仕様」の章に明記されているよう に、周囲温度がシェーカーの動作速度の範囲外になります。OKをタップして通知を確 認し、続行するか、別の温度を選択します。

注 温度設定値画面は、図 33の左下欄に示すように、校正温度と校正中に調整され た補正値における読み取り値を示します。温度校正の手順については、126ページの 「温度校正」。

 メイン画面に戻るには、表示される「設定が保存されました」確認ダイアロ グを了解します。

注 温度調節機能付シェーカーは、設定温度到達の際に、温度オーバーシュートと呼ばれ る特定の動作をします。温度オーバーシュートとは、チャンバー温度が、設定温度よりわ ずかに高い(または低い)温度にオーバードライブされた後に設定温度に近づくことを意 味します。設定温度に近づいている間、タッチスクリーンにはこのオーバーシュート動作 が表示されません。その代わりに、設定温度に達するまでチャンバー温度の上昇(または 下降)が表示されます。

シェーカーの開始と停止

シェーカーを開始するには、開始ボタンをタップします。
 開始ボタンが停止ボタンに変わります。



シェーカーを停止するには、停止ボタンをタップします。

3. 2. 2. ステータス

シェーカーの状態が良好な場合は、タッチスクリーンディスプレイの「ヘルスス テータスと統計」エリアに緑色のハートのアイコンが表示されます (図 26の⑦および図 27の®)。緑色のハートのアイコンをタップする と、ステータス画面が開きます。ステータス画面には、前回の動作時

の総振とう時間や総電源投入時間など、シェーカーの動作統計情報が表示されま す。温度調節機能付シェーカーには、さらに冷却や加熱にかかった時間が表示さ れます。

Status: Good		2	×	Status: Good			×
Total Run Time		Last Run Time Ses	sion	Total Run Time		Last Run Time Ses	sion
Hours Shaking	232hrs	Hours Shaking	60brs	Hours Shaking	232hrs	Hours Shaking	60hrs
				Hours Powered	332hrs	Hours Powered	61hrs
Hours Powered	332hrs	Hours Powered	61hrs	Hrs tempering	168hrs	Hrs tempering	61hrs
			(j)				(i)

図 35: シェーカーに関する統計:開放型シェーカー(左)と温度調節機能付シェーカー(右)

<u>アラート</u>

アラートが発生すると、タッチスクリーンディスプレイの現行画面の上部に黄色 のニュースティッカー型バーが表示されます。さらに、音響アラーム が鳴ります。メッセージを2回スクロールすると、黄色のニュースティ ッカー型のアラートバーが消えます。その後、黄色の三角形のみが表 示され、そのシェーカーにおいて1つ以上の アラートが存在することを示しま す。三角形アイコンは、有効なアラート件数が表示された円(周囲が白で青色の 円)とともに表示されます。「情報とヘルスステータス」エリア(図 26の⑦お よび図 27の⑧)にある三角形アイコンをタップすると、現在有効なすべてのアラ ートのリストを表示する画面が開きます。図 36に示すように、最新アラートは 詳細を確認できるように拡大表示されます。リストをスクロールし、閲覧するリ スト項目をタップすると拡大し、詳細が表示されます。



アラート項目の横にあるチェックボックスをタップすると、有効なアラートを選 択できます。了解ボタンをタップすると、リストからアラートの消去を試みま す。すべてのアラートが消去されると、ステータスアイコンがデフォルトの緑色 のハートに戻ります。

アラーム

アラームが発生すると、サンプルや装置への損傷を避けるためにシェーカーは直 ちに停止します。その際、必ずタッチスクリーンのアラームを確認し (___)
てから操作を続けてください。

アラートが発生すると、タッチスクリーン・ディスプレイの現行画面 の上部に赤色のバーが表示されます。「情報とヘルスステータス」エリア(図 26の⑦および図 27の⑧)に音の波の印で囲まれた赤いアラームベルが表示され ます。加えて、警報アラームが鳴り続けます。

赤いアラームバーの下にあるニュースティッカーには、現在のアラームの概要が スクロール表示で表示されます。右側に「スヌーズ」ボタンが表示され、アラー ムを一時的にミュートすることができます。スヌーズの間隔内にアラームを解除 しないと、警報アラームが再び鳴ります。スヌーズ間隔の時間は、69ページの 「アラームとアラート」のセクションで記述されている手順に従って、設定で選 択できます。

「情報とヘルスステータス」エリア(図 26の⑦および図 27の⑧)にあるベルア イコンをタップすると、図 37に示すように、現在有効なすべてのアラートのリ ストを表示する画面が開きます。リストをスクロールし、閲覧するリスト項目を タップすると拡大し、詳細が表示されます。



図 37: アラームのリスト

アラーム項目の横にあるチェックボックスをタップすると、有効なアラームを選 択できます。根本原因が解消されている場合は、了解ボタンをタップすると、リ ストからアラームの消去を試みます。すべてのアラームが消去されると、ステー タスアイコンがデフォルトの緑色のハートに戻ります。

<u>エラー</u>

誤動作が発生すると、シェーカーはエラーメッセージを表示し、サンプルやユニット自体の損傷を防ぐためにすぐに停止します。その場合、画面がすべて赤に変わり、以後操作不能になります。図38の例で示すように、エラーメッセージがエラーコードとともに表示されます。



図 38: エラー画面

問題を以下の手順で処理し、動作を再開してみてください。

- 1. エラー画面に表示されるエラーコードに注意してください。
- 2. 了解をタップして、アラーム音をミュートします。
- シェーカーの電源をオフにしてから再度オンにして、シェーカーを再起動し ます。
- それでもエラーメッセージが表示される場合は、Thermo Fisher Scientific カスタマーサービスに連絡し、エラー画面に表示されるエラーコードを提示 してください。

3. 2. 3. セッティング

ナビゲーションバー上の2番目の項目は、設定アイコンです。設定をタップする と、以下の画面が表示されます。設定画面には、表示ウィンドウに一度に表示で きるよりも多くのボタンがあります。図 39の右側に示すように、スクロールし て残りのボタンを表示してください。

<	Shake Lab 1 11:22:24AM Settings		<	Shake I 11:22:2 Setting	Lab 1 4AM s		
\$. €	Alarms and Alerts	î	¢, (8	Controls		Ŷ
Ë	U Access Control		Ë	Ģ	Display		
Ē	Calibration		Ē		Files and Info		
Ļ	a Controls	~	Ļ	×	Service	6	~
		0					()
図 39	設定画面						

アラームとアラート

「アラームとアラート」画面では、アラームとアラートの発行の時間と環境につ いて設定できます。



図 40: アラームとアラートにおけるプレファレンスの設定

音量: 下図の音量コントロールで、アラームやアラートが発生したときに鳴る音のスピーカーの音量が変更できます。スライダーを左右にドラッグし、音量を上 げ下げして調節し、保存をタップします。ここでの音量設定はアラームとアラー トの両方に適用されます。



図 41: アラームとアラートのスピーカー音量の設定

高温アラーム・低温アラーム(温度調節機能付シェーカーのみ): 設定温度に 対してそれぞれ高温・低温のアラームの閾値が設定できます。(64ページの「 温度設定」参照)。シェーカーのチャンバー内の温度が低温度閾値を下回わった 場合、または高温度閾値を超えた場合に、シェーカーがアラームを発行します。 温度閾値を選択し、保存をタップします。



注 設定温度を変更すると、それに応じてアラーム閾値が移行します。

高温シャットダウン・低温シャットダウン(温度調節機能付シェーカーのみ): 設定温度に対してそれぞれ高温・低温の安全停止の閾値が設定できます。(64 ページの「温度設定」参照)。シェーカーのチャンバー内の温度が低温度閾値を 下回わった場合、または高温度閾値を超えた場合に、シェーカーが自動的に運転 を停止して、チャンバーで処理中の試料を保護します。閾値を選択し、保存をタ ップします。

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

図 43: 高温・低温シャットダウン閾値の設定

注 設定温度を変更すると、エラー閾値が移行します。

スヌーズタイムアウト:スヌーズタイムアウトの時間では、メインウィンドウ上 部に表示される赤いバーのスヌーズボタンをタップしたときに、アラームが消音 される時間を決定します(67ページの「アラーム」参照)。スヌーズタイムア ウトの設定は5分、10分、15分から選択できます。デフォルトのオプションは10 分です。変更を保存するには、保存をタップします。

Snooze Timeout	5min ㅅ
	10min
	15min

図 44: アラームのスヌーズタイマーの設定

校正通知を無効にする(温度調節機能付シェーカーのみ):シェーカーの温度測 定を校正するように通知が定期的に表示されるのを無効に設定できます(126ペ ージの「温度校正」参照)。校正実行の通知を完全に無効にするか、プログラム 実行中に限り無効にするかを選択できます。

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

図 45: 校正通知を無効にする

アクセス制御

「アクセスコントロール」画面では、アクセス要件が制限されている手順に対応 するための管理者パスコードを割り当てることができます。

デフォルトでは、シェーカーは「オープンモード」になっています。この場合、 装置にアクセスして操作するためにパスコードを入力する必要はありません。「 セキュアモード」に切り替えると、シェーカーの設定を変更する際に管理者パス コードが必要になります。タッチスクリーンディスプレイの右上隅にログインボ タンが表示されます。

下記に、パスコードログインが必要な場合の「オープンモード」と「セキュアモ ード」の比較を示します。

アクション	オープンモードで必要 とされるパスコード	セキュアモードで必要 とされるパスコード
速度、時間、温度の設定値の設定	いいえ	いいえ
シェーカーの開始および停止	いいえ	いいえ
軌道の計算	いいえ	いいえ
アラームやアラートの了解	いいえ	いいえ
「ファイルと情報」画面の表示	いいえ	いいえ
ヘルスステータスの表示	いいえ	いいえ
イベントログとチャートの表示お よびエクスポート	いいえ	いいえ
アラームのスヌーズ	いいえ	いいえ
プログラムの実行	いいえ	いいえ
プログラムの作成、編集、削除	いいえ	はい
ディスプレイ設定の変更	いいえ	はい
コントロール設定の変更	いいえ	はい
アラームとアラートの設定の変更	いいえ	はい
アクセスコントロールの設定の変更	いいえ	はい
プログラムのインポートまたはエ クスポート	いいえ	はい
ファクトリーリセットの実行	いいえ	はい
校正の実行	いいえ	はい
ファームウェアのアップグレード のインストール	はい	はい

表 52: オープンモードおよびセキュアモードでのパスコードのログイン要件

図 46は、アクセスコントロール画面でのオープンモードとセキュアモードを示 しています。



図 46: アクセス制御: オープンモード(左)とセキュアモード(右)

セキュアをタップしてから保存をタップすると、変更を確認する前に、管理者パ スコードを入力するように表示されます。すべてのユニットは、工場で事前に設 定され、マニュアルに印刷されているものと同一の管理者パスコードとともに出 荷されています。

これと同じように、パスコードで保護された設定を変更する際には、管理者パス コードを入力するように表示されます。

注 事前に設定されている管理者パスコードは「00000」です。

管理者パスコードの変更ボタンで、事前に設定されたパスコードの変更ができま す。管理者パスコードの変更ボタンをタップして、最初に現在のパスコードを入 カし、次に新規パスコードを入力し、もう一度新規パスコードを入力して確認す るように表示されます。新規パスコードが保存されましたメッセージが表示され たら、変更手順が完了しています。

校正(較正)

「校正」画面(温度調節機能付シェーカーのみ)では、正確に温度を調節するために、シェーカーの校正を行う一連の画面を開くことができます。手順については、126ページの「温度校正」を参照してください。
コントロール

「コントロール」画面には以下のようなオプションがあり、ユニットで開始する デフォルトの動作パラメーターをあらかじめ設定することができます。

設定速度: 速度設定値を15rpmから525rpmの間に設定することができます。ホ イールピッカーのダイヤルを回転して速度を設定し、保存をタップします。「設 定速度」画面に関する使用方法の詳細な手順については、61ページの「設定速 度」のセクションを参照してください。

時間モード: シェーカーは連続モードまたはタイマーモードのどちらかで実行 できます。連続モードの場合は、ユーザーの判断でシェーカーを手動で停止しま す。タイマーモードの場合は、タイマーが切れるときにシェーカーの駆動が自動 的に停止します。タイマーモードでは、デフォルトのランタイム(運転時間)を 設定したり、シェーカー実行中のGUIの時間表示を経過時間か残り時間か選択し たりすることができます。「時間モード」画面に関する使用方法の詳細な手順に ついては、63ページの「ランタイム(運転時間)の設定」のセクションを参照 してください。

温度設定値(温度調節機能付シェーカーのみ): 起動時にシェーカーに表示す るデフォルトの温度を設定することができます。「温度設定値」画面に関する使 用方法の詳細な手順については、64ページの「温度設定」のセクションを参照 してください。

オートリスタート: この機能は、通常の開始・停止、プログラム、温度校正の実 行中に停電が発生した場合に回復後、ユニットを再起動します。オートリスター トがいいえに設定されている場合は、停電後、装置は自動的に再起動しません。

ディスプレイ

ディスプレイ設定を使用すると、さまざまな表示オプションが変更できます。 明るさ:ディスプレイの明るさレベルを調整するには、スライドコントロールま たは「+/-」ボタンを使用します。



図 47: 画面の明るさの調整

言語: 表示言語を変更するには、言語ボタンをタップします。ホイールピッカー を回転して使用する言語を選択し、保存をタップします。



図 48: 表示言語の選択

測定単位(温度調節機能付シェーカーのみ):測定単位をタップすると、GUI画 面全体の温度表示を°C(摂氏)と°F(華氏)の間で切り替わります。



図 49: 温度表示の単位の選択

日付:日付を設定するには、「日付」ボタンをタップします。ホイールピッカー の月、日、年のダイヤルを回転して選択し、保存をタップします。



図 50: 日付(月、日、年)の設定

日付フォーマット: 日付フォーマット を設定するには、「日付フォーマット」 ボタンをタップします。選択した日付形式に対応するラジオボタンをタップし (例:月/日/年(西暦)は4月15日2015年と表示されます)、保存をタップしま す。



図 51: 日付フォーマットの設定

時間: 時間と時間フォーマットを設定するには、「時間」ボタンをタップしま す。12時間または24時間をタップして、ホイールピッカーの時、分、12時間フォ ーマットの場合はAM/FMのダイヤルを回して、保存をタップします。



図 52: 時刻(12時間フォーマットまたは24時間フォーマット)の設定

スリープモード:スリープモードをタップして「オン」にすると、15分間操作し ないと、シェーカーのディスプレイがスリープ状態になります。スリープモード の場合は、画面にタップしてスリープ解除プロンプトが表示されます。

Off	
	Off

図 53: スリープモードを有効にする場合

地域:ユニットを使用する地域を特定するには、地域ボタンをタップします。「 都市、国」テキストボックスをタップして、都市名の最初の文字を入力します。 最初の3文字を入力すると、一致した提案が表示されます。一致するものを選択 するか、続けて完全な名前を入力して、オンスクリーンキーパッドで保存をタッ プします。



図 54: 地域を設定する

ユニット名: シェーカーに名前を付ける場合や、シェーカー名を変更する場合 は、ユニット名ボタンをタップします。ユニット名テキストボックスをタップし て、入力します。完了したら、オンスクリーンキーパッドで保存をタップしま す。



図 55: シェーカーに名前を付ける場合

メニューのカスタマイズ: メニューのカスタマイズボタンをタップして、メ インナビゲーションバーの下部にある2つのアイコンをカスタマイズします。目 的のアイコンを、メイン画面のエリアから交換するアイコンの上にドラッグしま す。次へをタップして確定します。 注 メニューのリセットをタップすると、ナビゲーションバーをいつでも工場出荷時の設 定に戻すことができます。



図 56: ナビゲーションバーのカスタマイズ

<u>ファイルと情報</u>

この画面には、シリアル番号、GUIアプリケーションソフトウェアとそのバージ ョン、シェーカーのメインコントローラーのファームウェアとそのバージョン、 パラメーターファイルとそのバージョン、現在のオペレーティングシステムが表 示されています。

ファクトリーリセットをタップして、シェーカーを工場出荷時のデフォルト設定 に戻すこともできます。ファクトリーリセットを実行するには、管理者パスコー ドが必要です。また、実行すると、グラフィカルユーザーインターフェイスを使 用して設定した全部の設定が消去されます。イベントログは、ファクトリーリセ ットを行っても削除されません。

< *	Shake Lab 1 11:22:24AM Files and Info	< *	Shake Lab 1 11:22:24AM Files and Info		
¢. (Serial Number 123TSX990960 H.M.I. 4.1.0 Main Controller 7000 0612 V3 Parameters 50154439 V1	¢, ∢ ≝	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	>	1 1
	0				()

図 57: ファイルと情報

さらに下スクロールすると、図 57の右側に示すように、ファームウェアボタン が表示されます。このボタンをタップすると、シェーカーの新しいファームウェ アをインストールできます。新しいファームウェアのインストールの詳細な手順 については、130ページの「4.7.ファームウェアのインストール」を参照して ください。

注 新しいファームウェアは、認定サービス技術者から入手してください。

アフターサービス

「サービス設定」は認定を受けたサービス技術者に限りアクセスすることができ ます。

3. 2. 4. プログラム

ナビゲーションバーでプログラムアイコンをタップすると、プログラムのリスト の表示や新規のプログラムの作成ができます。ここでは、プログラムの作成、編 集、削除、インポートおよびエクスポートを行うことができます。図 58は、新 品のデバイス(左)と以前から使用されているデバイス(右)の「プログラム」 ウィンドウを示しています。使用されているデバイスには、ユーザーが作成した プログラムのリストが表示されています。



図 58: プログラム

プログラムのリストの各項目には、プログラムの詳細を含む「プログラムクイッ クビュー」画面を開く目のアイコン ●と、プログラムの編集のために開く鉛筆の アイコン√があります。

全プログラムにおけるプレファレンスの設定

「プログラム」画面内で、詳細設定ボタンをクリックすると、詳細設定画面に移 動します。この画面では、プログラムを開始したときの初動操作が設定できま す。このプレファレンスは、作成するすべてのプログラムに適用されます。



図 59: プログラムにおけるプレファレンスの設定

1. プレファレンスボタンをタップします。

- タイマーの表示のオプションを選択して、以下のうち、プログラムのタイマ 一の表示方法を選択します。
 - > 経過時間:スタートボタンを押してからシェーカーが稼働している時間
 > 残り時間:タイマーが切れるまでシェーカーが動作し続ける時間
- プログラム開始時にシェーカーの実行をすぐに開始する場合は、プログラム を読み込んでシェーカーを開始オプションを有効にします。デフォルト動作 では、プログラムを実行するために開始ボタンをタップする必要がありま す。

注 プログラムを読み込んでシェーカーを開始動作は、シェーカーのフードが閉じ ている場合に限り機能します。フードを閉じるとすぐに、自動的にランが開始されま す。

4. 保存をタップします。

プログラムの作成

最大99個のプログラムを作成して保存できます。

- 1. 「新規作成」ボタンをタップします。
- 2. プログラム名を入力します。



図 60: プログラムの作成

 プログラムする温度(温度調節機能付シェーカーのみ)、速度、時間と分で ランタイム(運転時間)を入力します。
 温度調節機能付シェーカーの温度制御を無効にするには、温度フィールドを

空欄にしておきます。 4. プログラムにさらにステップを追加するには、スクロールダウンしてステッ プの追加ボタンをタップします。

5. 保存をタップしてプログラムを保存します。

プログラムの編集

- 1. 編集するプログラムの横にある鉛筆アイコンズをタップします。
- 必須フィールドを編集します。保存をタップします。プログラムは新たな変 更事項と共に保存されます。
- 画面下部のステップの追加を選択するとステップが追加できます。3つ以上のステップがある場合は、画面をスクロールしてステップの追加ボタンを表示してください。
 - 注 複数のステップを伴う振とう運転中に、アラームやエラーメッセージが発生した場合、シ ェーカーは自動的に停止します。アラートが発生した場合は、振とう運転は続行されます。



図 61: ステップの追加

プログラムの削除

- 編集するプログラムの横にある鉛筆アイコンをタップします。
- 2. 削除をタップします。削除の確認メッセージが表示されます。



図 62: プログラムの削除

3. 削除をタップします。選択したプログラムが削除されます。

プログラムの実行

- 1. 既存のプログラムを実行するには、実行するプログラムを選択します。
- プログラムのクイックビューを閲覧する場合は、プログラムの横にある目のアイコンのをタップします。
- 3. 「プログラムのクイックビュー」ウィンドウが表示されます。



図 63: プログラムクイックビュー

- ホーム画面でランをタップし、プログラムを読み込みます。
 選択したオートランの設定に従って、プログラムは直ちに開始するか、ホーム画面で開始ボタンをタップすると開始します。
- プログラムが完了すると、下図のウィンドウが表示されます。OKをタップします。



図 64: プログラム完了

プログラムのインポート

あるシェーカーで作成したプログラムを別のシェーカーにインポートすることが できます。これを行うには、まずシェーカーのUSBポートにUSBドライブを差し込 みます。

下の画面は、USBドライブで識別されたすべてのプログラムを示しています。



図 65: インポートするプログラムの選択

インポートするプログラムを選択します。インポートをタップします。

プログラムが正常にインポートされると、下図のウィンドウが表示されます。OK をタップします。ここでUSBは取り外すことができます。



図 66: プログラムのインポート完了

プログラムのエクスポート

シェーカー間でプログラムをエクスポートすることができます。その際、まず USBドライブが接続されていることを確認してください。

USBが接続されている場合は、下図の画面が表示されます。



図 67: エクスポートするプログラムの選択

エクスポートするプログラムを選択します。エクスポートをタップします。 エクスポートが完了すると、下図のウィンドウが表示されます。OKをタップしま す。ここでUSBは取り外すことができます。



図 68: プログラムのエクスポート完了

イベントログ

ナビゲーションパネルの3番目のタブはイベントログで、ユーザーイベントとシ ステムイベントの記録が保存されています。メニューバーで「イベントログ」ア イコンをタップすると、図 69に示すように、「イベントログ」画面が表示され ます。



図 69: イベントログ

この画面には、最近のイベントのリストとともに各イベントに関する日付とタイムスタンプが表示されます。

リストから項目を選択すると、個々のイベントの追加情報を表示することができます。

イベントは、下記のカテゴリに従って種類別にフィルターをかけることができます。

- アラーム
- アラート
- 設定の変更
- 開始/停止の操作
- プログラムの実行
- フードを開ける操作(温度調節機能付シェーカーのみ)

フィルターを選択すると、図 70に示すように、右側の「すべて表示」ボタンが 「フィルター・オン」に変わります。



図 70: フィルターをかけたイベントログ

イベントログのエクスポート

エクスポートするイベントドロップダウンリストから、エクスポートするイベントを選択します。ログ・レポートをエクスポートする形式を選択します。



図 71: エクスポートするイベントログの選択

日付範囲を所定の選択肢から選択するかカスタマイズします。

注 エクスポートできるのは、過去6か月のイベントに限られます。

 なお、エクスポートはUSBドライブでも実行できます。ログ・レポートを保 存するためのUSBドライブを挿入します。ログ・レポートをダウンロードす るには、ログのエクスポートボタンをタップします。



図 72: エクスポート用のUSBドライブを挿入

チャート

チャートには、時間経過に伴う速度や温度(温度調節機能付シェーカーのみ) に関するデータのグラフが表示されます。X軸は時間、Y軸は速度または温度を表 示します。

温度チャートを用いて、設定値、実際の温度、周囲温度を経過時間に対してプロ ットすることができます。編集ボタンでは、3本の曲線のどれを温度チャートに 含めるかを選択できます。さらに、表示範囲は、1日、7日、またはユーザー指定 の日数に変更することができ、各日については全24時間か選択した時間のいずれ かが選択できます。 リフレッシュボタンでは、設定に従ってグラフの表示が更新でき、チャートを最 後に読み込んでから、または更新してから記録した新しいデータが表示されま す。

注意:温度アラームが発生した場合、GUIに表示される温度は直ちにシェーカー チャンバーの実際の気温に切り替わり、サンプルを最大限に保護します。これに より、チャート機能に表示されるチャンバー温度に急激な変化が生じる場合があ ります。



図 73: 速度チャートおよび温度チャート

USBドライブをシェーカーのUSBポートにはめ込み、チャートデータのエクスポー トをタップすると、速度や温度のグラフのデータをダウンロードできます。

チャートの編集

1. 編集をタップして、チャートを編集します。



図 74: チャートの編集:記録する温度または速度の選択

- 温度または速度をタップして、表示するチャートを選択します。
- 3. さらに下スクロールして、日付範囲オプションを表示します。
- タップして、1日、7日、またはカスタム日付範囲の場合はカスタムを選択し ます。
- 5. ホイールピッカーのダイヤルを回して開始日を選択し、開始日を設定ボタン をタップします。

 ホイールピッカーのダイヤルを回して開始日を選択し、終了日を設定ボタン をタップします。

< *	Shake Lab 1 11:22:24AM Edit Chart			•	Custom Date Range From: Mar 14 2014 April 15 2015
¢.	Date Range:	1 Day7 Days		Î	May 16 2016
		From:	April 13, 2015		Custom Date Range To:
Ļ		lo: Sa	April 16, 2015 ve	•	April 15 2015 May 16 2016 Set From Date

図 75: チャート作成における日付範囲の選択

- 7. さらに下スクロールして、時間範囲オプションを表示します。
- タップして1日24時間を選択するか、カスタムをタップして特定の期間をチャートに記して、以下の操作をします。
 - a. ホイールピッカーのダイヤルを回して毎日の開始時刻を選択し、開始時 刻を設定ボタンをタップします。
 - b. ホイールピッカーのダイヤルを回して毎日の終了時刻を選択し、終了時 刻を設定ボタンをタップします。

注 開始時刻と終了時刻を選択することにより、2時点の間の正確な期間に対するチャートの作成を選択したことになります。1日あたり24時間のオプションを使用する 場合は、チャートは毎日00:00に開始し、23:59に終了します。



図 76: チャートを作成する時間帯の選択

9. 変更またはカスタム設定を保存するには保存をタップします。

3. 3. 付属品



注意

容器の破損による身体的・生物学的危害。 付属品が正しく取り付けられていないと、ガラスが割れたりサ ンプルが流出したりすることがあります。 付属品が正しい工具とネジで適切に取り付けられているか確認 してください。 付属品がプラットフォームに適切に組み合わさっているか確認 してください。

容器は必ず適切な寸法の付属品で使用してください。



注意

鋭利な端部による切り傷。

プラットフォームなどの付属品の取り扱いには注意してくださ い。



注意

プラットフォームの動作中に指がはさまれる恐れがあります。 動作中は、絶対に指をプラットフォームの上下に置かないでく ださい。

注 ユーザーは責任を持って、付属品が正しく設置されているか確 認してください。

取り付け方法は、本取扱説明書に別途記述がない限り、すべてのシェーカーモデ ルに適用されます。

常に正しい工具を使用し、特に付属品については供給されているものを使用して ください。付属品を紛失したときのために、スペアパーツキットを注文すること ができます(17ページの「付属品」)。指定されている工具以外のもの、また供 給されているネジ以外のものは使用しないでください。

3. 3. 1. プラットフォームの設置

注意 各シェーカーに対応するプラットフォームの完全なリストについては、18ページ の「1.2.1.プラットフォーム(振とう台)」セクションを参照してください。 温度調 節機能付シェーカーモデルについては、工場で事前に取り付けられたプラットフォームに 限り動作できます。

注意 鋭利な端部による切り傷。プラットフォームを取り外すときは、プラットフォーム の下側をつかみます。プラットフォームに取り付けてあるクランプでプラットフォームを 持ち上げないでください。プラットフォームなどの付属品の取り扱いには注意してください。

注意 Solaris 4000 I/4000 Rシェーカーで11x14プラットフォームを動作させると、手 や指が挟まり、押しつぶされることがあります。Solaris 4000 I/4000 Rシェーカーでは 11x14プラットフォームを使用しないでください。

注意 プラットフォームには付属のネジとロックツール以外を使用しないでください。他 のネジや不適切な固定工具を使用すると、取り付け不備の原因になり、シェーカーや付属 品が損傷する可能性があります。

必ずすべてのネジを使用してプラットフォームを取り付けてください。 必ず適切なロックツールを使用してください。

- 3/16インチロックツール (ネジ締め) (GT530066) : Solaris 2000、 Solaris 2000 I/Rおよび4000 I/R シェーカー用
- 7/32インチロックツール (ネジ締め) (GT530080) : Solaris 4000シェー カーの全プラットフォーム 用

<u>標準プラットフォーム</u>

Solaris 2000および4000の各シェーカーには、標準プラットフォーム1台、ネジ 数個、ロックツール1つが付属しています。アプリケーションに合わせて追加の プラットフォームが購入できます。各シェーカーに対応しているプラットフォー ムの完全なリストについては、18ページの「1.2.1.プラットフォーム(振 とう台)」セクションを参照してください。

2000 I、2000 R、4000 I、4000 Rシェーカーには、シェーカーにプラットフォー ムがあらかじめ組み込まれています。プラットフォームやネジなどの個々の部品 はありません。



図 78: シェーカーに取り付けられた標準プラットフォーム: Solaris 2000 (左)、Solaris 4000 (右)

 プラットフォームを慎重に取り付ける位置に合わせて、シェーカーの上に水 平に置きます。

Solaris2000シェーカーのプラットフォームには、取付穴が3つありま す。Solaris 4000シェーカーのプラットフォームには、取付穴が4つあります。 Solaris 2000 I/Rおよび4000 I/Rシェーカーのプラットフォームには、取付 穴が3つあります。

ネジをゆっくりと締めて、プラットフォームをシェーカーに固定します。ロックツールが曲がり始めたら、締めるのをやめます。

<u>デュアルスタック・プラットフォーム</u>

注 デュアルスタックプラットフォームは、Solaris 2000 I/Rおよび4000 I/Rシェーカー には対応していません。



図 79: デュアルスタック・プラットフォームの組立て

上段プラットフォームと下段プラットフォームを各隅にある4本のピラーを使用 して組立てます。上段プラットフォームと下段プラットフォームを適切なネジを 使用してピラーで取り付けます。

ネジをゆっくりと締めて、ピラーとプラットフォームを接続します。ロックツー ルが曲がり始めたら、締めるのをやめます。

シェーカー接続用の取付穴がある方が下段プラットフォームです。組み立てると きは、必ず下段プラットフォームを下側にしてください。 注 必ず組立てたプラットフォームがぐらつかないか最終チェックを行ってから、その上 に容器を置くようにしてください。



図 80: Solaris 4000シェーカーのデュアルスタック・プラットフォーム

- デュアルスタック・プラットフォームを慎重にプラットフォーム取付板に合わせて、シェーカーの上に水平に置きます。
 Solaris 2000用のプラットフォームには3つの取付穴があり、Solaris 4000 用のプラットフォームには4つの取付穴があります。
- ネジをゆっくりと締めて、プラットフォームをシェーカーに固定します。ロ ックツールが曲がり始めたら、締めるのをやめます。

3. 3. 2. フラスコクランプと容器の取付け

注意 容器の破損・液漏れによる生物学的危害。付属品が正しく取り付けられていないと サンプルが流出する可能性があります。付属品が正しい工具とネジで適切に取り付けられ ているか確認してください。取付けには、シャフトの長さが15cm(6インチ)のドライバ ー (PH2)が必要です。(商品番号 75004131)。付属品がプラットフォームに適切に組み合 わさっているか確認してください。容器は必ず適切な寸法の付属品で使用してください。 容器は損傷がないものを正しく取り付けてください。

注意 鋭利な端部による切り傷。プラットフォームなどの付属品の取り扱いには注意して ください。

注意 クランプを安全に固定するには、付属のネジとの組み合わせが必要不可欠です。付 属のネジを、そのクランプの指定された取付穴全部に取り付けます。

注 ネジが長すぎると、交換可能なプラットフォームの安定性と機能が損なわれる可能性 があります。固定するには、付属品に含まれているネジのみを使用してください。

92 / 138

<u>クランプ</u>

各クランプには、容器クリップ1個、スプリング1個または2個(クランプ・ボト ルのサイズによる)、プラットフォームに取り付けるためのネジ(複数)が含ま れています。付属のネジ以外はクランプに使用しないでください。



スプリングチューブ ② スプリング ③ ネジ ④ クリップレッグ ⑤ スプリング
 図 81: 2個のスプリングとクランプの詳細

- 1. 必要に応じて、図のようにスプリングをクリップレッグに取り付けます。
- 図に示されているように、ゴム製のスプリングチューブがクリップレッグの 間になるようにセットします。クランプには2つのスプリングを使用するも のもあります。2つ目のスプリングをプラットフォームに取り付けてから、 それを組立てたクリップの基部周辺に取り付けます。
- 3. 組立てたクリップを付属のネジでプラットフォームに取り付けます。

<u>容器</u>

- 使用する容器をクランプに配置する際、注意して最初に クランプのスプリングを十分に引き離して容器の基部が クランプの内側に配置します。容器をゆっくりと適切な 位置にスライドさせて、クランプのより広い底部に固定 します。スプリングで容器の首をしっかりと固定し、所 定の位置に保持します。
- すべての容器がしっかりと固定されているか確認して、 装置の電源を入れます。

混合動作中に物質が散乱しないように、可能な限り、容 器をストッパーで密閉する必要があります。



3. 3. 3. 角型クランプの取付け

注意 容器の破損・液漏れによる生物学的危害。付属品が正しく取り付けられていないと サンプルが流出する可能性があります。付属品が正しい工具とネジで適切に取り付けられ ているか確認してください。取付けには、シャフトの長さが15cm(6インチ)のドライバー (PH2)が必要です。(商品番号 75004131)。付属品がプラットフォームに適切に組み合わ さっているか確認してください。容器は必ず適切な寸法の付属品で使用してください。容 器は損傷がないものを正しく取り付けてください。

注意 鋭利な端部による切り傷。プラットフォームなどの付属品の取り扱いには注意して ください。

94 / 138

<u>角型クランプの詳細</u>

角型クランプは、パネル底に取付穴とあり、ばね仕掛けのクリップレッグ4つが 付いた構造になっています。また、プラットフォームにクランプを取り付けるた めのネジも含まれています。付属のネジ以外はクランプに使用しないでくださ い。



① 取付穴 ② スプリング ③ クリップレッグ ④ ネジ

図 82: 角型クランプの詳細

- 必要に応じて、図 82に示すように、すべてのスプリング②をクリップレッグ ③に引っ掛けます。
- 2. クリップアセンブリをプラットフォームに配置し、取付穴を合わせます①。
- 3. 付属のネジ④で、クリップアセンブリをプラットフォームに取り付けます。
- 4. 上から四角い容器を挿入します。
- 5. 容器が底部のパネルに接触していて、クランプにしっかりと固定されている ことを確認します。

3. 3. 4. 試験管立ての取付け

注意 鋭利な端部による切り傷。プラットフォームなどの付属品の取り扱いには注意して ください。

試験管立ては、固定ノブ付取付ブラケット(ワイヤーフレームまたはシートメタ ル)2個と、プラットフォームに取り付けるためのつまみネジ2つが各試験管台に 付いた構造になっています。試験管立てアセンブリには、試験管立てに発泡イン サートがあらかじめ組み込まれています。必ずラックに付属のネジを使用してく ださい。



① 試験管立て
 ② 発泡体インサート
 ③ つまみネジ
 ④ 取付ブラケット
 ⑤ 固定ノブ

図 83: ワイヤーフレーム取付ブラケット付きの試験管立てアセンブリ



図 84: シートメタル取付ブラケット付きの試験管立てアセンブリ

- つまみネジで、取付ブラケットをプラットフォームに取り付けます。手でつ まみネジをできるだけ締めます。
- 2. 試験管立てを取付ブラケットにはめ込みます。
- 固定ノブ⑤で必要な角度に設定します。手で固定ノブをできるだけ締めます。

3. 3. 5. 角度調節式試験管立てホルダーの取付け

注意 鋭利な端部による切り傷。プラットフォームなどの付属品の取り扱いには注意して ください。

試験管立てホルダーと試験管立ては、28ページの「1.2.5.角度調節式試験 管立てホルダー」に記載されているような、さまざまな寸法のものが利用可能で す。すべての試験管立てホルダーは7つの位置に調整可能で、0°から両方向にス イングし、15°、30°、45°でロックします。

必ずラックホルダーに付属のネジを使用してください。



図 85: 試験管立てを取り付けた角度調節式試験管立てホルダー

- 付属のネジで、ラックホルダー①をシェーカーのプラットフォームに取り付けます。
- ラックホルダーのスイングベッド⑥の両端にある金属製のタブ③を少し広 げ、試験管立て②をはめ込みます。
- 3. 慎重にタブ③を放して、試験管立て②をしっかりと固定します。
- ホルダーの両端にあるロッキングピンのノブ⑤を外側に引きます。ノブを4 分の1回転すると、ピンのロックが解除されます。
- ラックホルダー①のスイングベッド⑥を15°、30°、45°のうち使用する角度に回転させます。
- ロックピン⑤を位置決め穴④にもう一度取り付け、ノブを4分の1回転してピンをロックします。
- ラックを取り外すには、スイングベッド⑥の両端にある金属製のタブ③を広 げ、試験管立て①を持ち上げます。

 3. 6. マイクロプレート・ディープウェルプレートのクラン プの取付け

注意 鋭利な端部による切り傷。プラットフォームなどの付属品の取り扱いには注意して ください。



ディープウェルプレート
 ネジ
 マイクロプレートクランプ

図 86: マイクロプレート・ディープウェルプレートのクランプアセンブリ

- 1. マイクロプレートのフレームをプラットフォームに置きます。
- マイクロプレート・ディープウェルプレートセットに付属のネジで、マイク ロプレートのフレームをプラットフォームに取り付けます。
- マイクロプレートまたはディープウェルプレートをマイクロプレートのフレ ームに挿入します。
- マイクロプレートやディープウェルプレートを静かに持ち上げて、しっかり と取り付けられていることを確認します。

3. 3. 7. 調節式容器クランプの取付け

注意 鋭利な端部による切り傷。プラットフォームなどの付属品の取り扱いには注意して ください。





① 取付穴 ② 可動ロッド ③ 固定ノブ ④ トレイの側壁 ⑤ 滑り止めマット ⑥ ネジ

図 87: 調節式一段式容器クランプアセンブリ



6 7 7 7 7

① 取付穴 ② 可動ロッド ③ 固定ノブ ④ トレイの側壁 ⑤ 滑り止めマット ⑥ ネジ

図 88: 調節式二段式容器クランプアセンブリ

以下の手順で取り付けます。

- 滑り止めマットを持ち上げて、取付穴(図 88に示す①)が見えるようにします。
- 調節式容器クランプをプラットフォームに配置し、取付穴を位置合わせます。
- クランプに付属のネジ⑥で、調節式容器クランプをプラットフォームに取り 付けます。
- 4. 滑り止めマット⑤を調節式容器クランプの底に敷きます。
- 5. 容器を滑り止めマット⑤の上に置きます。
- 6. 固定ノブ③を緩めます。
- 可動ロッド②の位置を調節して、容器をしっかりと固定します。
- 8. 固定ノブ③を締め直します。
- 9. 容器を静かに持ち上げて、しっかりと固定されていることを確認します。

100 / 138

3. 3. 8. ビーカーラックの取付け

注意 鋭利な端部による切り傷。プラットフォームなどの付属品の取り扱いには注意して ください。



図 89: ビーカーラックアセンブリ

- 1. パネル底②をプラットフォームの上に載せます。
- 2. ビーカーラック①をパネル底の上に載せます。
- 3. 取付穴をプラットフォームに合わせます。
- 4. 製品に付属のネジ③で、ビーカーラックアセンブリを固定します。

3. 3. 9. 分液ロートクランプの取付け

注意 鋭利な端部による切り傷。プラットフォームなどの付属品の取り扱いには注意して ください。

注 必ず正しいネジを使用してください。付属のネジは2種類あります。赤いスレッドロッ カー付きの長い方のネジは、垂直スタンドの固定に使用します。スレッドロッカーのない 短い方のネジは、分液ロートクランプをプラットフォームに取り付ける場合に限り使用し ます。

分液ロートクランプを用いて、分液ロートを基本的に2方向、つまり水平および 垂直に、シェーカーのプラットフォームに取り付けることができます。水平およ び垂直。



図 90: 分液ロートクランプの取付位置

図 90の水平位置③で高さを調整すると、④に示すように傾斜した向きにすることができます。さらに、分液ロートクランプを垂直スタンド(付属品として供給されている)と組み合わせると、①および②に示すように垂直に配置することができます。また、クランプは分液ロートの栓を①のように下に向けて取り付けるか、反転して②のように逆さまに配置することができます。

垂直取付の準備

図 91は、分液ロートクランプのアセンブリとその付属品の垂直スタンドを横並 @びに示してい®ます。



 ④ 分液ロートクランプ
 ⑧ 垂直スタンド
 ① プラットフォーム取付用 平らな穴
 ② ロートクランプ取付用ネジ筋付穴
 ③ 標準垂直スタンド用取付穴
 ④ 反転垂直スタンド用取付穴
 ⑤ 赤いスレッドロッカー付きネジ
 ⑥ スレッドロッカーなしの ネジ

図 91: 垂直スタンド付属品(右)付分液ロートクランプ(左)

- 1. 図 90の①および②に示すように、分液ロートクランプを垂直の取付台の長い 方の端に取り付けます。
- 2. 垂直スタンドのロートクランプ取付穴4つ(図 91の②)に、以下の手順で位 置合わせします。
 - a. 標準垂直取付穴(図 91の③)を使用して、分液ロートの栓を下に向けて 取り付けます。
 - b. 反転垂直取付穴(図 91の④)を使用して、分液ロートの栓を下に向けて 取り付けます。
 - c. 分液ロートクランプを付属のネジ(図 91の⑤)で、垂直取付台に固定 します。

プラットフォームへの取付け

- 1. 分液ロートクランプアセンブリをプラットフォームに配置します。
- 2. 取付穴をプラットフォームに合わせます。
- 分液ロートクランプに付属のネジ(図 91および図 92の⑥)で、分液ロート クランプアセンブリをプラットフォームに固定します。

103 / 138

<u>分液ロートの取付け</u>



① 分液ロート
 ② 後方ロートホルダー
 ③ 前方ロートホルダー
 ④ 高さ調整用固定ノブ
 ⑤ 長さ調整用固定ノブ
 ⑥ スレッドロッカーなしのネジ

6

図 92: 分液ロートクランプアセンブリ

- 1. 長さ調整用固定ノブ⑤を緩め、前方ロートホルダー③を前方の端までスライ ドさせます。
- 2. 分液ロート①の後端を後方ロートホルダー②にはめ込みます。
- 前方ロートホルダー③を後方にスライドし、分液ロートの脚を前方ロートホ ルダー③の開口部に配置します。
- 前方ロートホルダー③を適切な高さに調整し、高さ調整用固定ノブ④を締めて固定します。高さを調整することにより、分液ロートを傾斜した位置に傾けることができます。
- 5. 長さ調整用固定ノブ⑤を締めて、前方ロートホルダー③を固定します。

3. 3. 10. ユーティリティトレイの取付け

注意 鋭利な端部による切り傷。プラットフォームなどの付属品の取り扱いには注意して ください。



図 93: ユーティリティトレイアセンブリ

以下の手順で取り付けます。

- 1. ユーティリティトレイ②をプラットフォームの上に載せます。
- 2. 取付穴をプラットフォームに合わせます。
- ユーティリティトレイに付属のネジ③で、ユーティリティトレイ②を固定し ます。
- 4. ゴム製マット①をトレイ②に敷きます。
- 5. 図 94に示すように、振とうするボトルをトレイに置き、付属の0リング④で 固定します。
 - 注 100rpmを超える速度で振とうする場合は、0リングでの固定が推奨されます。



図 94: 培養容器が0リングで取り付けられたユーティリティトレイ

105 / 138

注 容器を固定するときは、0リングを伸ばしすぎないように注意してください。0リ ングを取り付けるために伸ばすと、そこを超えると0リングが伸びなくなる点、つま り弾性における静止点があることに気が付きます。この静止点に達したら、0リング が切れるため、それ以上伸ばさないでください。

3. 3. 11. 粘着マットの取付けと使用

粘着マットは底が広く平らな容器、つまり三角フラスコ、フェルンバッハフラス コ、培地ボトル、マイクロプレート、シャーレー、細胞培養フラスコ、メスフラ スコ、ビーカーなどに適しています。一般的に直立可能なすべての容器を使用で きますが、底面積と高さの比が低下すると、達成可能な最大の振とう速度も低下 します。図 95に示すように、底が小さく高い容器はあまり適していません。大 きな三角フラスコ(例:3000ml)は、小さなフラスコ(例:100ml)よりも強く 付着します。



図 95: 粘着マット: 適切なフラスコの形状

粘着マットについては、以下の2つの一般的な種類が利用可能です。

- 低接着マット(透明) 品番75004111、75004117
 - » 粘着性が少なく扱いやすい
 - » ビーカーに最適

» 最高200rpmまで(組織培養による増殖用)のすべての容器に適している

- 高粘着マット(黒色透明) 品番75004126、75004127
 - » 粘着性が高く容器を確実に固定する
 - » すべての容器に適している
 - » 最高速度400rpm (500ml までの三角フラスコ使用)

道具類

図 96に示す道具は、ローリングアプリケーターと呼び、粘着マットの付属品として供給されています。これは、粘着マットをプラットフォームに押さえつけるために使用します。



図 96: 粘着マット用ローリングアプリケーター

ドライマットの取付け

- 1. シェーカーのプラットフォームが清潔で乾燥していることを確認します。
- 2. 粘着マットの片側から保護ホイルを剥がします。
- 剥がした粘着面を下にしてマットをプラットフォームにしっかりと貼り付け ます。
- 4. 上面の保護ホイルを取り外します。

ウェットマットの取付け

- 1. 粘着マットの両面から保護ホイルを剥がします。
- 2. マットを水に浸すか、水でよくすすぎます。
- 3. マットから水滴が落ちるのを待ちます。
- 湿ったマットを清潔なプラットフォームに広げ、必要に応じて位置を合わせ ます。
- プラットフォーム上でマットを24時間乾燥させてから使用します。

容器の特性に関する留意事項

最高速度は容器の種類によって異なり、容器の大きさ、形状、底の形状、素材、 表面粗さ、清浄度などによって分類されます。シェーカーを使用する前に、これ らのパラメーターを評価して、用途に最適なマットを選択する必要があります。 表 53を用いて、粘着マットの使用に関する素材別実験器具の適性を判断しま す。

プラスチック/ガラス器具	粘着度	金属	粘着度
ガラス(ホウケイ酸ガラス)	5	ステンレス鋼	5
ポリプロピレン (PP)	3	アルミニウム	5
ポリプロピレン共重合体 (PPC0)	3	陽極酸化アルミ ニウム	4
ポリエチレン(HPDE) (LPDE)	3		
ポリカーボネート (PC)	4		
ポリエチレンテレフタレート (PET)	4		
ポリエチレンテレフタレート コポリエステル(PETG)	4		
ポリメチルペンテン (PMP)	3		
ポリテトラフルオロエチレン (PTFE)	0*		
テトラフルオロエチレン・ヘ キサフルオロプロピレン共重 合体(FEP)	0*		
パーフルオロアルコキシポリ マー (PFA)	2		
シリコーン	1*		
評価: 0 = まったく粘着性がな	い、5=最た	大の粘着性を示す/ * 評	価が0および1の容

器(PTFE、FEP、シリコン)は使用しないでください。

表 53: 素材別容器の適性

注 最高振とう速度は、表 53で「5」と評価された素材に限り達成可能です。

注 ガラス以外の素材については、素材別に実際の負荷、速度、ランタイム(運転時間)、 温度で実行できるかを各自で試して決定する必要があります。
付属品

ガラス容器(30%充填時)における速度範囲および運転時間に対する制限 ガラス容器(30%充填時)における室温での速度および運転時間に対する制限。 温度を上げると粘着性が減少します。温度が変更する手順では、容器がマットから離れる可能性があります。容器の温度が順化した後、振とうを開始します。

	マイクロプレート・ ディープウェルプ レート	ビーカー			
	96 x 2 ml	25 ml	125 ml	250 ml	500 ml
125 rpm	運転時間の制限なし				
150 rpm	72時間	72時間	72時間	72時間	72時間
200 rpm	72時間	72時間	72時間	72時間	72時間
250 rpm	72時間				
300 rpm					
350 rpm					
400 rpm					

表 54: 低粘着マット - 室温での速度範囲と運転時間に対する制限

	マイクロプレー ト・ディープウ ェルプレート	エルレ	ンマイヤ	ー フラ	スコ				
	96 x 2 ml	25 ml	125 ml	250 ml	500 ml	1000 ml	2000 ml	3000 ml	5000 ml
125 rpm	運転時間の制限なし								
150 rpm	72時間	72時間	72時間	72時間	72時間	72時間	72時間	72時間	72時間
200 rpm	72時間	72時間	72時間	72時間	72時間	72時間	72時間	72時間	72時間
250 rpm	72時間	72時間	72時間	72時間	72時間	72時間	72時間	48時間	24時間
300 rpm	72時間	72時間	72時間	72時間	72時間	72時間	48時間	24時間	
350 rpm	48時間	48時間	48時間	48時間	48時間	48時間	24時間		
400 rpm	24時間	24時間	24時間	24時間	24時間				

表 55: 高粘着マット - 室温での速度範囲と運転時間に対する制限

準備

注意 ガラス器具を取り扱うときは、必ず保護メガネと手袋を使用してください。 粘着マットに容器を乗せる前に、以下の手順を行います。

- 1. 容器に損傷がないか確認し、損傷しているものは使用しないでください。
- 清潔で乾いたワイプで容器を拭いて、水分を完全に取り除きます。
- グリースなどの汚染物質を完全に取り除きます。
- 4. 容器の底と粘着マットが清潔で乾燥した状態であることを確認します。

注意 どんなに少量の水であっても、容器の底に残っていると容器との接着を緩め、 しばらく動作した後、マットから外れる可能性があります。



図 97: 粘着マット: 容器を拭いて乾燥する

5. 液体の流出を防ぐために 、充填量を下げるか、容器を密閉してください。

注意 容器が完全に付着している場合でも、一滴の液体が外壁をつたって流れ落ちる と、しばらく動作した後、マットから外れる可能性があります。

容器の取付け

- 1. 容器をマットの上に置き、押し下げます。
- 軽く押したり引いたりすると、容器がマットに確実に接着します。

注 低温または長時間の運転では、凝結水が形成して、容器が離れる事故が発生す る可能性がありますのでご注意ください。試料と容器の温度は、シェーカーの指定 温度範囲から逸脱してはいけません。シェーカーの可能な温度範囲の詳細について は、11ページの「1.1. 仕様」を参照してください。 容器の取外し

- ディスプレイウィンドウにある「停止」ボタンを押すか、ランが終了するの を待ちます。
- プラットフォームが完全に停止するまで待ちます。
- 容器の首の部分をゆっくりと均等に横に引き、数秒待ちます。容器の底面が 大きい場合は、取外しに数秒かかる場合があります。



図 98: 粘着マット:容器を慎重に押したり引いたりします

注意 物理的な力をあまりかけずに、軽く押したり引いたりします。容器が突然緩ん で、内容物が流出する可能性があります。

4. 容器が外れない場合は、容器の周囲にごく少量の水を垂らします。

再びマットを使用する前に、風乾します。

粘着マットの取外し

 粘着マットを外すには、図 99に示すように、片側を持ち上げてプラットフ ォームからゆっくりと引き離します。



図 99: 粘着マット:マットを持ち上げて引き剥がして、取り外します

<u>クリーニングと消毒</u>

粘着性が低下したら、マットを清掃してください。

注意 スキンケア添加剤を含まない中性食器用洗剤のみを使用し、それ以外の溶剤、洗 剤、洗浄剤は使用しないでください。洗剤の推奨濃度は、水1リットルあたり最大2mlで す。 注 濃縮洗剤がマットに直接接触すると、マットの粘着性が減少します。その場合、すす ぎだけでは不十分です。洗剤が直接掛かってしまった場合は、マットをすすぎ、24時間水 に浸してから、24時間乾燥させて完全な接着力を回復します。

注 70%イソプロパノールまたは70%エタノールのみで消毒します。

わずかに汚れている、またはほこりが付いている場合

- プラットフォーム上のマットを上側からウェットワイプで拭くか、消毒します。
- 2. マットを湿らせます。
- 3. スポンジで残留物を拭き取ります。
- 4. 浄水に浸したスポンジで水拭きします。
- 5. 必要に応じて消毒します。
- 6. 24時間乾燥させます。

ひどく汚れている場合

- 完全に汚れを落とすために、プラットフォームからマットを取り外します。
- 希釈した食器用洗剤入りの水にマットを浸します。
- スポンジでマットを隅々までこすり、すべての残留物を取り除きます。
- きれいになったら、マットを水で十分に洗い流します。
- 必要に応じて消毒します。
- 6. 24時間乾燥させます。
- 107ページの「ドライマットの取付け」または107ページの「ウェットマットの取付け」の説明に従って取り付けます。

3. 3. 12. ガス供給マニホールドの取付け

ガス供給マニホールドを使用すると、シェーカーの使用中に窒素や二酸化炭素な どの不活性ガスを複数の容器に注入できます(注文する場合は、42ページの 「1.2.12.その他の付属品」を参照してください)。

ガス供給マニホールドは、Solaris 2000 I/Rおよび4000 I/R シェーカーには出 荷時に取り付けられていますが、それがないユニットには、ユーザー自身で取付 可能な付属品としてもお求めになれます。Solaris 2000 I/Rおよび4000 I/Rに はどちらかのサイドパネルにアクセスポートがあり、そこから実験用ガス供給チ ューブでマニホールドアセンブリにガスを供給できます(ポートの位置について は、44ページの「1. 4. 製品概要」を参照してください)。ガス供給マニホ ールドには、ガスを供給する入口継手1つと、容器に対する出口継手8つがありま す。



図 100: Solaris 4000 R シェーカーに取り付けられたガス供給マニホールド

<u>以下の手順で取り付けます。</u>

- 付属のネジ⑤で、ガス供給マニホールド②をシェーカーチャンバーの後壁に しっかりと固定します。
- アクセスポートのハトメの外側からプラスチックのプラグを取り外します (位置については、44ページの「1.4.製品概要」を参照してください)。
- シェーカーチャンバー内にあるゴム製ハトメの膜にカッターナイフで、交差
 印(X)の切り込みを入れます。
- 切断可能なチューブをカットして、実験用ガス供給チューブ①として使用するのに適した長さにます。

- 5. 実験用ガス供給チューブ①をマニホールドの逆目付きの入口継手④に取り付 けます。
- 実験用ガス供給チューブ①がフードのすべてのレバーから離れていることを 確認し、アクセスポートのゴム製のハトメを通して外側に送ります。
- 7. 実験用ガス供給チューブ①をガス供給源に取り付けます。
- 8本のチューブを切り取り、マニホールドの出口③に取り付けます。
 容器が9個以上ある場合は、Yコネクターを使用します。容器が7個以下の場合は、ガスの損失を削減するために使用しないチューブは挟んで止めます。
- 9. 適切なプラグとシールを使用して、チューブを容器に送ります。

3. 4. 負荷と 通常操作



化学反応の誘発による火災の危険性。

シェーカーは容器の中身が散乱するような速度で運転しないで ください。

警告 速度はゆっくり上げてください。化学薬品を使用するときは、 まず水で試してください。



装填の不備や付属品の損傷は安全性を低下させる恐れがあります。

注意

- 付属品とサンプルを含めた負荷がプラットフォームの中心に 対して対称に配置されているか確認してください。デュアル スタック・プラットフォームの操作時に、負荷重量が上段プ ラットフォームよりも下段プラットフォームの方が重くなっ ているか確認してください。
- シェーカーに過大な負荷をかけないでください。最大負荷については、(11ページの「仕様」)を参照してください。
 負荷重量には、シェーカーに取り付けられているプラットフォーム、付属品、およびサンプルの重量が含まれます。
- シェーカーを操作する前に、付属品が正しく取り付けられているか確認してください。88ページの「付属品」の指示に従ってください。



フードが落ちて閉じると、指や手が負傷する危険があります。

注意

シェーカーのフードが不意に閉まって、指や手が挟まった り、圧迫されたりして、負傷することがあります。これは、 フードのスプリングに欠陥があり、フードが完全に開いた位 置で保持できない場合に発生することがあります。 フードのスプリングに欠陥がある状態でシェーカーは使用し ないでください。Thermo Fisher Scientificのカスタマー サービスに連絡して、フードのスプリングをすぐに交換して ください。 読み込む

付属品やサンプルを含む負荷がプラットフォームの中心に対して対称にかかるように配置されているか確認してください。デュアルスタック・プラットフォーム を運転する際に全容量を使用していない場合は、 最高の性能を実現するため に、できれば底板に負荷をかけてプラットフォームの中心に対して対称的に負荷 がかかるように配置します。

温度調節機能付シェーカーを使用するときは、積載物の損傷や衝突を防ぐため に、フードを閉じるときに十分な間隔を確保し、フードを閉じた状態で適切に振 とうできるように、試料が置かれていることを確認してください。



図 101: 負荷が適切なプラットフォームの例



図 102: 負荷が適切でないプラットフォームの例

注 最大負荷重量には、シェーカーに取り付けられているプラットフォーム、付属品、お よびサンプルの重量が含まれます。最大負荷の詳細については、11ページの「1.1.仕 様」を参照してください。 通常操作

速度負荷の制限に関する情報は、製品の使用上の手引きを目的としてのみ提供さ れています。低振動または無振動での円滑な運転は、充填量、クランプやマット の種類と状態、容器の種類、プラットフォームの容器の配列、シェーカーの設置 表面、設定速度などの様々な要因によって異なります。

新規または未知の組み合わせを設定するときやこれらのパラメーターのいずれか を変更するときは必ず、 速度をゆっくりと175rpm以上に上げ、シェーカーの負 荷に対する不要な振動や動作を確認してください。ユーザーは責任を持って、シ ェーカーを安全に操作してください。

Solarisシェーカーは、特に低熱出力かつ低振動に重点を置いて設計されている ため、環境室、インキュベーター、実験用冷蔵庫などのさまざまな実験機器内で の使用に非常に適しています。実験機器内で使用する場合は、棚や固定補助具に 関する安定性など、さまざまな特定の環境条件が異なるため、安全な操作に対す る責任はユーザー自身が負うものとします。ゆっくりと速度を早めて、不要な機 器の動作を確認します。

安全なスピード

Solarisシェーカーに対する負荷非依存の安全速度については、表 56を参照して ください。負荷重量には、シェーカーに取り付けられているプラットフォーム、 付属品、およびサンプルの重量が含まれます。最大負荷の詳細については、11 ページの「仕様」を参照してください。

モデル	最高安全速度
Solaris 2000	175 rpm
Solaris 4000	175 rpm
Solaris 2000 I	250 rpm
Solaris 2000 R	250 rpm
Solaris 4000 I	250 rpm
Solaris 4000 R	250 rpm

表 56: 最高安全速度

最高速度

Solarisシェーカーでは、最高速度525rpmで作動できる負荷が表 57のように限られています。負荷重量には、シェーカーに取り付けられているプラットフォーム、付属品、およびサンプルの重量が含まれます。使用するクランプと容器が高速度には適さないこともあるので、それらの速度性能を考慮に入れてください。

モデル	最大安全負荷
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

表 57: 最高速度での最大負荷

注 速度が上がるとクランプが開き始め、容器の動きが大きくなるため装置に一時的な振 動が加わります。

重量対速度曲線

下記の重量対速度の曲線は、通常の使用に最適な速度と負荷の組み合わせを設定 するときのガイドです。緑色で示された領域は、振とう運転中に振動が発生しな い速度と負荷の組み合わせを示しています。赤色で示された領域は、振とう運転 中に強い振動が発生し、シェーカーの不要な動きを招く可能性がある速度と負荷 の組み合わせを示しています。これらのチャートはあくまでも目安ですので、速 度と負荷の組み合わせの設定が赤色の領域に接近している場合は注意する必要が あります。

ユーザー独自のアプリケーション条件によっては、表示された境界線に到達する 前に不要な負荷や装置の動きが生じることがありますので注意してください。負 荷に対する動き具合を調べるために速度は徐々に上げてください。負荷重量に は、シェーカーに取り付けられているプラットフォーム、付属品、およびサンプ ルの重量が含まれます。

注 以下の各グラフで青色の水平線は、機種別のプラットフォームの空虚重量を示してい ます。



図 103: Solaris 2000 - 通常操作



Solaris 4000

図 104: Solaris 4000 - 通常操作



図 105: Solaris 2000 I / 2000 R - 通常操作



Solaris 4000 I / 4000 R

図 106: Solaris 4000 I / 4000 R - 通常操作

4. メンテナンスとケア



警告

有害物質の取り扱いによる危険

- 有害物質を振とうさせる際は、世界保健機関(WH0)の 「実験室バイオセーフティマニュアル」および現地の規 定を遵守してください。世界保健機関(WH0)の「実験 室バイオセーフティマニュアル」に従って、リスク群2 の微生物サンプルを振とうさせる場合は、エアロゾル封 じ込め用バイオロジカルシールを必ず使用して用バイオ ロジカルシールを必ず使用してください。世界保健機関 WH0のホームページ(www.who.int)にて、「実験室バイ オセーフティ指針」をご検索ください。価の高いリスク グループの物質の場合、予防対策を必ず二つ以上備えて ご利用ください。
 - シェーカーやシェーカーの部品が毒素や病原物質に汚染 された場合は、適切な消毒対策を取る必要があります (124ページの「除染」、124ページの「消毒」の参 照)。
 - 危険な状況が発生した場合は、シェーカーの電源を切り、速やかにその場を離れてください。



警告

感染性物質による健康被害

事故による流出などで液体や物がプラットフォームの下に入 り込んだ場合は、直ちにシェーカーの電源を切り、プラグを 抜き、プラットフォームを取り外してください(18ページ の「プラットフォーム(振とう台)」)。通常の検査法に従 って、流出物を清掃してください。適切な個人用保護具を使 用してください。

装置内の調整や修理は、必ず認定を受けたサービス技術者に要請してください。 ユーザーはシェーカーのハウジングを開けてはいけません。

適切な保管および清掃方法については、付属品に付いてくる製品情報に記載され ている具体的な詳細に従ってください。本章の以下の情報はガイドラインとして のみ使用してください。

4. 1. 基本事項

個人保護、環境保護、および機材保護のため、シェーカーと付属品は必ず定期的 に清掃し、必要に応じて消毒してください。

Thermo Fisher Scientificは、お使いのシェーカーは最低でも毎月1回清掃し、 手作業で消毒することを推奨します。標準的な室内の空気中には何千もの微生物 が循環しており、それらが、シェーカー内に留まることによって、培養が危険な 状態にさらされることがあります。

Thermo Fisher Scientificは、70%エタノール、70%イソプロパノール、または 10%以下の第四級アンモニウム塩系の消毒剤の使用を推奨します。

注 振動機構に液体が入ると故障することがあります。液体、特に有機溶剤は振動機構や ベアリングに接触させないようにしてください。有機溶剤に触れると、ベアリングのグリ ースが取り除かれます。クリーニングできないシェーカーの部分に液体が染み込んだ場合 は、シェーカーを動作せず、フィールドサービスに連絡してください。

注 評価されていない手順や化学物質を用いると、シェーカーの素材を劣化させ、誤動作 や故障の原因になる可能性があります。意図した手順が機器に安全であるという確信がな い限りは、その他の洗浄や汚染除去の手順の実行は避けてください。機器に損傷を与えな い洗浄剤のみを使用してください。不確かな場合は、クリーニング剤のメーカーにお問い 合わせください。

注 温度調節機能付シェーカーでは、洗浄剤と水が蒸発するように、フードを開けたまま にしておきます。

- まず、電源プラグを抜き、クリーニング、消毒、除染を行います。
- シェーカーに取り付けられている付属品とプラットフォームを取り外してから、クリーニング、消毒、除染を行います。
- 中性洗剤を温水に溶かし、柔らかい布に含ませて拭き清掃します。ご質問が あれば、Thermo Fisher Scientificまでお問い合わせください。清浄水で洗 い流した後、完全に乾燥させます。
- リン酸、漂白剤、クレンザーなどの腐食性洗浄剤は絶対に使用しないでください。
- 消毒剤はpH値が6~8のものを使用します。
- 液体が流出した場合は直ちに、非腐食性の洗浄剤で湿らせた毛羽立ちのない 布を使用して、洗浄剤の製造元の指示に従って清掃します。
- 流出液はプラットフォームの下に染み込むことがあります。プラットフォームの下にこぼれたら、プラットフォームを取り外し、プラットフォームの下にあるトイレから流出液を完全に取り除きます。必要に応じて、消毒または除染します。
- シェーカーの部品を確認し、流出した培養液や残屑を取り除きます。

付属品の点検

注 破損の兆しのあるシェーカーや付属品は使用しないでください。安全性を確保するた めに、日常点検の一環として、定期的な付属品の点検を推奨します。

完全に付属品を清掃したら、損傷や摩耗、腐食がないか検査してください。

金属部品

腐食、磨耗、亀裂などの損傷が見られた付属品は直ちに使用を中止してくださ い。

プラスチック部品

ひび割れ、退色、傷、亀裂などがないかを確認してください。検査で破損が見ら れた製品は、直ちに使用を中止してください。

粘着マット

粘着マットを使用する前に必ず、摩耗や損傷の兆候がないか検査します。たとえ ば、ひび割れ、退色、黄変、もろさ、変形、表面の擦り傷、薬品による腐食など を確認します。粘着マットに摩耗や損傷の兆候が見られた場合は、すぐに廃棄し ます。

4. 2. クリーニング

注 クリーニングを実行する前に、実行予定の方法が機器に損傷を与えないか洗浄剤メー カーに問い合わせて確認してください。

注 振動機構に液体が入ると故障することがあります。振動機構やベアリングに液体、特 に有機溶剤を付着させないでください。有機溶剤に触れると、ベアリングのグリースが取 り除かれます。 中性洗剤と水の混合液で、ユニットの外側を柔らかい布で洗い、清浄水で 洗い流し、完全に乾かします。

シェーカーと使用後の付属品の適切なクリーニングについては、122ページの「 基本事項」を参照してください。

注 温度調節機能付シェーカーでは、洗浄剤と水が蒸発するように、フードを開けたまま にしておきます。

タッチスクリーン

タッチスクリーンのクリーニング:

- 1. 電源コンセントを引き抜きます。
- タッチスクリーンは、乾いたマイクロファイバーの布を使用してきれいにし ます。
- 必要ならば、水で湿らせたマイクロファイバーの布で、もう一度タッチスク リーンを拭いてください。
- 4. 風乾させます。

123 / 138

4.3. 消毒

警告 有害物質の取り扱いによる危険。汚染された部品には触れないでください。汚染されたパーツへの接触は感染の危険性があります。容器が割れたり液が流出したりして、感染性物質がシェーカー内に入り込むことがあります。汚染が発生した場合は、誰も危険にさらされていないことを確認してください。該当パーツの消毒を直ちに行ってください。

注 不適切な消毒方法や薬剤の使用によって機器が破損することがあります。消毒剤や手 順が機器に損傷を与えないことを確認してください。確信が持てない時は、消毒剤の製造 元にお問い合わせください。ご使用の消毒剤についてはその安全上の注意事項および取り 扱い説明書に従ってください。

- 必要な消毒レベルに従って、70%エタノールですべての部品と領域を拭きます。暴露した電子機器の部分は濡らさないでください。
- 1. 風乾させます。

注 温度調節機能付シェーカーでは、洗浄剤と水が蒸発するように、フードを開けた ままにしておきます。

責任を持って、適用される要件に従って消毒レベルを達成してください。

4. 4. 除染

警告 有害物質の取り扱いによる危険。汚染された部品には触れないでください。汚染された部品に接触すると、暴露する危険性があります。チューブが割れたり液が流出したりして、汚染物質がシェーカー内に入り込むことがあります。汚染が発生した場合は、誰も危険にさらされていないことを確認してください。該当のパーツをすぐに除染してください。

注 不適切な除染方法やクリーニング剤による、機器の破損。除染剤や手順が機器に損傷 を与えないことを確認してください。ご質問があれば、除染剤の製造元にお問い合わせく ださい。ご使用の除染剤についてはその安全上の注意事項および取り扱い説明書に従って ください。

Thermo Fisher Scientificは以下の方法を提案します。

- 70%エタノールですべての部品と領域を拭きます。暴露した電子機器の部分 は濡らさないでください。
- 2. 風乾させます。

注 温度調節機能付シェーカーでは、洗浄剤と水が蒸発するように、フードを開けた ままにしておきます。

責任を持って、適用される要件に従って除染レベルを達成してください。

4. 5. オートクレーブ

注 オートクレーブ温度と時間に関する許容値を、絶対に超えないようにしてください。 蒸気での化学品の添加は、許容されていません。

高圧蒸気滅菌処理の準備:

すべての付属品を分解します。

化学薬品や洗剤をよく洗い流します。最後に蒸留水ですすぎます。
 高圧蒸気滅菌処理が可能か次の表で確認してください。

品目	オートクレー ブ可・不可	仕様
シェーカー	いいえ	-
プラットフォーム	はい	121°C、20分
クランプ(スプリングチューブなし)	はい	121°C、20分
ねじ回し	いいえ	-
マイクロプレート・ディープウェルプレートクランプ	はい	121°C、20分
粘着マット	いいえ	-
ビーカーラック	はい	121°C、20分
分液ロートクランプ	はい	121°C、20分
試験管立て - 金属部品	はい	121°C、20分
試験管立て - プラスチック部品	いいえ	-
試験管立て - 発泡インサート	いいえ	-
調節式容器クランプ - 金属部品	はい	121°C、20分
調節式容器クランプ - 滑り止めマット	いいえ	_
ユーティリティトレイ - 金属トレイ	はい	121°C、20分
ユーティリティトレイ - ゴム製マット	いいえ	-
角度調節式試験管立てホルダー - 金属部品	はい	121°C、20分
角度調節式試験管立てホルダー - プラスチック部品	いいえ	-
角度調節式試験管立てホルダー - 発泡インサート	いいえ	-

表 58: 高圧蒸気滅菌処理の可能な素材

地域や施設などの要件に準じて、必要とされる無菌状態が達成されることを確認 してください。

高圧蒸気滅菌処理後、外観や色がわずかに変化することがあります。

4. 6. 温度校正

温度調節機能付シェーカーの設定温度は、精度が既知である温度測定装置に対し て校正して、恒温(加熱)または冷却(冷却・加熱)の振とう処理が完全に再現 できるようにすることができます。

4. 6. 1. 温度校正の最良実践法

シェーカーの温度校正を実行する前と実行中は、正確に手順に従っていることを 確認してください。推奨される最良実践法から外れたり逸脱したりすると、サン プル温度が不安定になり、校正の実行が失敗する可能性があります。

最良実践法は以下の手順で行います。

- 250mlの容量のフラスコを使用します。
- フラスコに正確に100mlの水を入れます。
- フードを完全に閉じ、校正を実行している間は常に、閉じたままにします。
- +分な時間をとって、サンプル温度が完全に安定するまで待ちます。
- 校正の実行中に、温度や湿度などの環境条件が変化しないようにします。ドアや窓の開閉や空調システムの設定の変更など、小さなことでも、継続している温度校正の手順の結果に支障を来す可能性があります。
- シェーカーの校正は年に1回行うことが推奨されます。

4. 6. 2. 温度の校正の手順

校正するには、下記の手順で行います。

- 1. 校正を開始するには、以下の3通りがあります。
 - » 54ページの「初回起動」に記載されている初回起動の手順の最後に、 初回起動の手順を終了する代わりに、温度校正を実行するオプションが 選択できます。
 - » 年に一回校正するようにシェーカーに通知が表示されます。
 - » メニューバーから設定アイコンをタップし、続いて表示される温度校 正画面のさらに下にある校正ボタンをタップすると、校正が開始できま す。
 - » メニューバーから設定アイコンをタップし、続いて校正ボタンをタップ すると、必要に応じていつでも手順が開始できます。
 - 注 校正の通知は、設定画面の校正通知を無効にするチェックボックスで無効にする ことができます。

図 107に示す温度校正画面が表示されるのを待ちます。



図 107: 校正:開始画面

3. 校正を開始するボタンをタップして開始します。

Shaker Lab 1 01:47:52PM Temp calibration
Step 1 of 5
Fill a 250-mL Erlenmeyer flask with 100mL of water and position it (fixed with a clip) at the
approximate geometric center of the shaking platform.
Next

図 108: 校正:三角フラスコに水を充填する手順の表示

- 画面に表示される手順に従います。250mlの三角フラスコに100mlの水を入れ ます。
- フラスコをクランプに取り付けて、このクランプをプラットフォームのほぼ 中央に固定します。
- 精度が既知である温度測定装置のプローブをフラスコに入れ、水中に浮かせ ます。プローブは水中に沈めますが、フラスコの底や側面に接触してはいけ ません。
- 7. 水がこぼれないようにフラスコを密閉します。
- 8. プラットフォームの上のフードを閉じます。

9. 次へボタンをタップして続行します。



図 109: 校正:校正を開始する手順の表示

注 開始ボタンをタップすると、シェーカーがすぐに振とうし始めます。

画面に表示される手順に従います。開始ボタンをタップして、あらかじめ設定した温度を使用するか、設定温度フィールドをタップして目的の温度を入力し、次に開始ボタンをタップします。

画面には、シェーカーのチャンバー内において、選択した校正温度への加熱 (または冷却)状況が表示されます。



図 110: 校正:チャンバーは校正温度に調整中

 チャンバーの温度測定が安定状態になるまで、最低でも2時間半(150分間) 待ちます。

次の図 111に示すように、安定状態になったら、チャンバー温度が安定しま したというメッセージが表示されます。



図 111: 校正:チャンバーは校正温度で安定しました

- 12. 画面に表示される手順に従います。
 - » 図 111に示す画面が表示されたらすぐに、手続きに従って、熱電対温度 計の数値を比較します。温度が安定していると思われる場合は、編集ボ タンをタップして修正するか、終了ボタンをタップして温度校正の手順 を完了します。



図 112: 校正:シェーカーが表示した正しい温度

- 表示されるホイールピッカーで、シェーカーが表示する温度を正確な測定装置の読み取り値に修正します。
- 14. 保存ボタンをタップして、変更を登録します。

15. 温度校正が完了しましたのメッセージは、手順が正常に完了したことを示し ます。



図 113: 校正:温度校正が完了しました

再校正をタップすると、温度校正の手順を繰り返すことができます。

4. 6. 3. 温度校正の探索

現行の校正温度はいつでも調べることができます。

- メニューバーから設定アイコンをタップし、続いて表示される温度較正画面 のさらに下にある校正ボタンをタップします。
- 図 107に示す温度校正画面が表示されます。この画面には、現在のシェーカ 一校正温度が表示されます。

注 また、温度調節機能付シェーカーの場合は、温度を設定した画面で校正温度を 調べることができます。64ページの「温度設定」のセクションに例が示されていま す。

4. 7. ファームウェアのインストール

「ファイルと情報」画面(77ページの「ファイルと情報」参照)で「ファーム ウェア」ボタンをタップすると、管理者パスコードの入力するように表示され、 次に、図 114に示す「ファームウェア」画面が開きます。この画面では、USBメ モリから新しいファームウェアをインストールできます。

サービス契約によっては、現場のサービス技術員から新しいファームウェアをダ ウンロードするリンクの通知があります。ファームウェアは正規の供給元からの みインストールしてください。承認されていない供給元からのソフトウェアを使 用すると、T Thermofisher Scientificの認めた保証が無効になる場合がありま す。

- 現場のサービス技術員から供給されたリンクからZIPファイルをダウンロー ドします。
- 2. ZIPファイルを空のUSBドライブに展開します。

130 / 138

注 USBドライブに、ZIPファイルを展開するのに十分なスペースがあることを確認してください。

- シェーカーが運転中の場合は、表示ウィンドウの停止ボタンを押して停止し ます。
- 66ページの「ステータス」のセクションで記述されている手順に従って、 現存するすべてのアラームを了解します。

注 シェーカーの運転中にアップグレードしようとすると、「シェーカーが振とうし ている間は、シェーカーのアップグレードができません」というメッセージが表示さ れます。そのメッセージには、シェーカーの停止ボタンがあり、運転を終了してアッ プグレードが実行できるようにします。

- 5. ナビゲーションバーのファイルと情報アイコンをタップします。
- ファイルと情報画面を一番下までスクロールして、ファームウェアボタンを 表示します。
- ファームウェアボタンをタップして、現在インストールされているファーム ウェアバージョンが表示されているファームウェア画面を開きます。
- USBドライブをシェーカーのUSBポートの1つに差し込みます。USBポートの位置については、44ページの「製品概要」のセクションを参照してください。
- 9. 管理者パスコードを入力します。



図 114: ファームウェア画面

10. ファームウェアのアップグレードボタンをタップします。

注 シェーカーがUSBドライブを認識できない場合は、「インストールするファーム ウェアを含むUSBメモリを差し込んで、 アップグレードを開始してください」という メッセージが表示されます。ここでUSBドライブを差し込むか、すでにUSBドライブが はめ込まれている場合は、USBポートに正しく取り付けていることを確認し、問題を 修正してください。 ファームウェアのインストール

 アップグレードするファームウェアがある場合は、図 115に示すような項目 別リストが「ファームウェアのアップグレード」画面に表示されます。アッ プグレードするファームウェアの項目をタップします。

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

図 115: インストールするファームウェアの項目の選択

注 最新のファームウェアの項目はグレー表示され、選択できません。更新する項目がま ったくない場合は、「現時点ではソフトウェアの新しいアップグレードはありません」と いうメッセージが表示されます。

- 12. 選択をタップします。
- USBメモリにファームウェアのバージョンが複数含まれている場合は、1つ選 択するように表示されます。ファームウェアのバージョンをタップして選択 し、選択をタップします。



図 116: インストールするファームウェアのバージョンの選択

注 どのバージョンをインストールするかがわからない場合は、ダウンロードするリンク を提供した現場のサービス技術者にお問い合わせください。 ファームウェアのインストール

アップグレードする項目のリストが表示されます。OKをタップしてアップグレードを開始します。



図 117: アップグレードする項目のリスト

注 GUIではソフトウェアの相互従属性を処理することができます。つまり、更新するファ ームウェアの項目を1つ選択すると、そのソフトウェアが必要とする全項目を自動的にイン ストールに追加します。

注 GUIファームウェアのみをアップグレードするため、図 115に示されているリストの オプションでH.M.I.を選択した場合は、アップグレードを開始してから最高1分間は、GUI ディスプレイでタッチスクリーンのアクティブなエリアをそれ以上タップしても応答しま せん。この間は、シェーカーの電源を<u>切らない</u>でください。電源を切った場合は、アップ グレードが正常に完了しません。シェーカーはしばらくすると再起動し、また使用できる ようになります。

注 複数のファームウェアのコンポーネントを一度にアップグレードするため、図 115に 示すリストから2つまたは3つのオプションを選択した場合は、アップグレードの処理中は 常にシェーカーの電源をオンに保持します。電源を切ってはいけません。

 図 118に示すように、インストールの進行状況を通知するさまざまなメッセ ージが表示されます。



図 118: ファームウェアをアップグレードするときのインストール中のメッセージ

シェーカーを再起動するように指示が表示されたら、電源スイッチでユニットの電源をオフにして、再びオンにします。

4. 8. プラットフォームの交換

プラットフォームの取付け方に関する基本的な手順は、89ページの「3.3.1. プラットフォームの設置」を参照してください。各シェーカーに対する予備のプ ラットフォームの完全なリストについては、18ページの「1.2.1.プラット フォーム(振とう台)」セクションを参照してください。

注 温度調節機能付シェーカーモデルについては、工場で事前に取り付けられたプラット フォームに限り動作できます。

4.9. アフターサービス

Thermo Fisher Scientificは、毎年一回、認定サービス技術者にシェーカーと 付属品の点検を要請することを推奨します。カスタマーエンジニアは以下の点検 を行います:

- 電子装置
- 使用場所の適切性
- 安全システム
- 使用した付属品
- クランプ、プラットフォーム、付属品などのシェーカーへの固定

サービスの予定日までに、シェーカーと付属品が完全に清掃、除染され、検査が 完全かつ安全に実行できる状態にしておいてください。

これらのサービスは、Thermo Fisher Scientificサービス・点検契約でご提供いたします。修理が必要な場合は、保証内であれば無料で、保証外の場合は有料で行います。これは、Thermo Fisher Scientificの認定サービス技術者がシェーカーのメンテナンスをしている場合に限り有効となります。

4. 10. 送付と廃棄

警告 感染性物質による健康被害。シェーカーと付属品を廃棄する場合は清掃し、必要に 応じて消毒・除染を行ってください。不明点がある場合は、Thermo Fisher Scientific カ スタマーサービスにご連絡ください。

シェーカーを廃棄する場合は各国の規定に従ってください。シェーカーの廃棄に 関するお問い合わせは、Thermo Fisher Scientificカスタマーサービスまでご連 絡ください。お問い合わせ先は本説明書の最終ページか、次のホームページをご 覧くださいwww.thermofisher.com。

運搬および運送についての記載もご覧ください(52ページの「運搬」ペ59ペ ージの「運送」参照)。 送付と廃棄

5. トラブルシューティング

数	説明	解決方法	アイコン
エラー			
11999	内部エラー発生。	内部エラーが発生しました。電源スイッチをオフにし てから再度オンにして、ユニットの電源を入れ直して ください。エラーメッセージの表示が消えない場合 は、サービスマンにご連絡ください。	
アラート			
3	温度校正の期日。 1年がたちました。	1年間の有効期間が終了しました。 126ページの「4.6.温度校正」のセクションで記 述されている手順に従って、温度校正を更新して ください。	CAL
4	電源が再投入され、オート リスタートが有効になり ました。	前回の実行中に停電が発生しました。電力の回復 後、実行が自動的に再開されました。	Ţ
6	周囲温度センサーが故障し ました。	周囲温度センサーによって測定された温度が不自 然です。	III
10	ファン1が最高回転速度を超 えています。	ファン1の回転速度が妥当速度の閾値を超えてい ます。	2
11	ファン2が最高回転速度を超 えています。	ファン2の回転速度が妥当速度の閾値を超えてい ます。	
12	ファン3が最高回転速度を超 えています。	ファン3の回転速度が妥当速度の閾値を超えてい ます。	2
13	ファン4が最高回転速度を超 えています。	ファン4の回転速度が妥当速度の閾値を超えてい ます。	
20	フードが長時間開いてい ます。	チャンパー温度を維持するためにシェーカーのフー ドを閉じてください。	
アラーム			
1	高温アラーム	チャンバー温度が最高アラーム値を超えています。 試料、環境条件を確認するか、設定を変更してく ださい。	
2	低温アラーム	チャンバー温度が最低アラーム値を超えています。 試料、環境条件を確認するか、設定を変更してく ださい。	
3	停電後のオートリスタート ができません。	前回の実行中に停電が発生しました。オートリスタ ートが正常に実行されません。	Ğ
5	ドライブの加速が遅すぎま す。 目的の設定速度に到達 できません。	時間内には目的の設定速度に到達できません。設定 やプラットフォームの負荷(付属品とサンプル)を 確認してください。	\bigcirc

135 / 138

数	説明	解決方法	アイコン
7	ファン1の速度計測が予期せ ぬ停止を示しています。	ユニットの電源を一旦切り、また入れてください。 エラーメッセージが続く場合は、サービス技術員に 連絡してください。	•2
8	ファン2の速度計測が予期せ ぬ停止を示しています。	ユニットの電源を一旦切り、また入れてください。 エラーメッセージが続く場合は、サービス技術員に 連絡してください。	2
9	ファン3の速度計測が予期せ ぬ停止を示しています。	ユニットの電源を一旦切り、また入れてください。 エラーメッセージが続く場合は、サービス技術員に 連絡してください。	2
10	ファン4の速度計測が予期せ ぬ停止を示しています。	ユニットの電源を一旦切り、また入れてください。 エラーメッセージが続く場合は、サービス技術員に 連絡してください。	•
12	ドライブの起動エラー発生 - 速度信号受信不可能。	シェーカーのブラットフォームがブロックされました。プラットフォーム周辺に十分なスペースがあるか確認してください。また、プラットフォームの負荷(付属品とサンプル)を軽減してください。その後、シェーカーの「開始」ボタンをもう一度押してください。エラーメッセージが続く場合は、サービス技術員に連絡してください。	٢
24	ラン処理中の速度計測エラ ーの発生。	異常な速度変化が検出されました。シェーカーのプ ラットフォームの負荷(付属品とサンプル)やクラ ンプを確認してください。その後、シェーカーの「 開始」ボタンをもう一度押してください。エラーメ ッセージが続く場合は、サービス技術員に連絡して ください。	
26	実行中に速度計測の予期せ ぬ停止が検出されました。	プラットフォーム周辺に十分なスペースがあるか確 認してください。また、プラットフォームの負荷(付属品とサンプル)を軽減してください。 シェーカーの背面にあるヒューズが切れていない ことを確認してください(47ページの「ヒュー ズ」参照)。 その後、シェーカーの「開始」ボタンをもう一度押 してください。エラーメッセージが続く場合は、サ ービス技術員に連絡してください。	٢
82	モーター電流測定のタイ ムアウトエラーが発生し ました。	モーターの過電流が検出されました。実行中にプラ ットフォームに負荷を加えたり減らしたりしないで ください。速度を落とすか、プラットフォームの負 荷 (付属品とサンプル)を調整してください。	— 0
83	モーターの電流が測定範囲 外です。	モーターの過電流が検出されました。実行中にプラ ットフォームに負荷を加えたり減らしたりしないで ください。速度を落とすか、プラットフォームの負 荷(付属品とサンプル)を調整してください。	— 0

表 59: エラー、アラート、アラームのリスト

注 この表にリストされていないエラーメッセージが表示された場合は、サービス技術員 に連絡してください。

GPL (General Public License:一般公有使 用許諾)

本機のソフトウェアの一部は、GPL、LGPL、または別のオープンソースライセン スに従って、公開されているオープンソースソフトウェアを使用しています。詳 しくは、これらは表にリストされているライブラリです。関連するライセンスに よって提供される場合、使用するライブラリ(サードパーティ製ライブラリ)の ソースコードはThermo Fisher Scientificから入手できます。使用するオープン ソースソフトウェアのそれぞれのライセンス条項は、提供されるソースコードパ ッケージの一部です。

サードパーティ製ライブラリ _____

ライブラリ	バージョン	展開する手段	ライセンス
Qt	5.8	BSP	LGPLv3
Log4Cplus	1. 2. 0	アプリ	Apache ライセン ス v2/2条項BSDラ イセンス
boost	1. 72. 0	アプリ	Boostライセンス 1.0
json11	1.0.0	アプリ	MITライセンス

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

thermo scientific

専用仕様





Thermo Electron LED GmbH Zweigniederlassung Osterode Am Kalkberg, 37520 Osterode am Harz Germany 原産国: USA

		Thermo	Scientific	Solaris	2000	Thermo	Scientific	Solaris	4000
(H	Thermo	Scientific	Solaris	2000 I	Thermo	Scientific	Solaris	4000 1
		Thermo	Scientific	Solaris	2000 R	Thermo	Scientific	Solaris	4000 F

70900190 は取扱説明書の原書です。 この取扱説明書は、オリジナルの取扱説明書を翻訳したものです。

thermofisher.com/shaker

© 2019-2024 Thermo Fisher Scientific Inc. 全権利を留保します。

すべての商標は、特記のない限り、Thermo Fisher Scientific Inc.およびその子会社の所有物です。国によっては、お取り扱いのない製品もございます。詳細は、お近くの代理店にお問い合わせください。

マニュアルに掲載されている写真は一例を示すものであり、パラメータや言語により異なる場合があります。

スペイン/ポルトガル +34 93 223 09 18

スイス +41 44 454 12 12

イギリス/アイルランド +44 870 609 9203

アメリカ合衆国・カナダ +1 866 984 3766

その他のアジア諸国 +852 2885 4613

その他の国々 +49 6184 90 6000

