thermoscientific



Thermo Scientific

STERI-CYCLE i160 LK / STERI-CYCLE i250 LK

CO2 インキュベータ

使用説明書

50162768 A

10 2 0 2 1

[©]2021 Thermo Fisher Scientific Inc. 全権利を留保します

商標

HERACELL VIOS™, Steri-run™, Steri-cycle™, iCan™, THRIVE™ および Cell locker™ は、Thermo Scientificの商標です。 Thermo Scientificは、Thermo Fisher Scientific Inc. 社の商標です。 本使用説明書において記載のある商標は、該当製造者の所有物です。

Thermo Electron LED GmbH Robert-Bosch-Straße 1 D - 63505 Langenselbold ドイツ

Thermo Electron LED GmbH社は、以下の子会社です: Thermo Fisher Scientific Inc. 168 3rd Avenue Waltham, MA 02451 USA

Thermo Fisher Scientific Inc.社は、本装置をご購入いただいたお客様に対し、装置使用のため 本書をご提供いたします。本書は、著作権で保護されています。Thermo Fisher Scientific Inc. 社の書面における承諾がない限り、たとえ一部であっても複写は禁止されています。

本使用説明書の内容は、予告なく変更される場合があります。本書に記載のある技術仕様は全て 情報のためであり、それに拘束力はありません。本書に記載のあるシステムコンフィギュレー ションおよび技術仕様は、場合によっては購入者が過去に得た情報を置き換えるものです。

Thermo Fisher Scientific社は、本書の完全性、正確性及び非欠陥性に対する請求は提起せす、 本書を使用した結果により生じ得るエラー、脱落、事後損傷については、例えこれらが本書記載 のデータに依るものであるとしても、その責任は一切負いません。

本書は、Thermo Fisher Scientific Inc.社と顧客との間で結ばれる購入契約書には含まれません。本書には、一般購入規約の内容に影響を及ぼすことはありません。一般購入規約と相違がある場合、一般購入規約の内容が優先となります。

目次

前書去	0–1
	0 1
一般事項	0-1
· 装置の情報及び記録	0-1
	0-1
利用者への指示	0-1
説明書の有効性	0-2
翻訳用の原本	0-2
保証	0-2
保証条件	0-2
安全上のご注意とマークの説明	0–3
本使用説明書における安全上のご注意とマークの説明	0–3
安全上の注意を示すマーク:	0–4
装置に記載のあるマーク	0-5
装置の利用目的	0–5
規定の利用目的	0–5
利用目的に反した使用	0-6
規定と要綱	0-6
ガスに対する安全上の注意	0-7
職員への指示:	0-7
二酸化炭素における安全上の注意(CO2)	0-7
酸素における安全上の注意(02)	0-8
窒素における安全上の注意 (N2)	0-8
装置	1–1
女世記之	1_1
已夜	1_1
11.20mmの11.11.11.11.11.11.11.11.11.11.11.11.11.	1-2
STERI CVCI E i160 LK / i250 LK 追加装備の配送内存	1_2
STEDI CVCI E 1160 LK / 1230 LK 迫加表開の配と内谷	12
SIENI-UIULIIOULN 追加表開の能区内谷	1-3
STERI-CYCLE 1250 LK 追加装備の配送内谷	1-3
装置の設置	2-1
周辺条件	2-1
西龙车佰·	2_1
—————————————————————————————————————	2 1
次 ス	2 Z 2_2
必安はへ、 へ	2 Z 2_2
建区 · · · · · · · · · · · · · · · · · · ·	2-3
波 l で 生 ね る	2-3
サホートノフケットの取り付け	2-0
里ね直さワアリエーション	2-1
可動式台座の取り付け	2-8
可動式台座と足(i160 LK と i250 LK での例)	2–9
追加装備、装備変更	2-13

1

2

3	装置の説明	3-1
	STERI-CYCLE i160 LK / i250 LK 前方から	3–2
	STERI-CYCLE i160 LK / i250 LK 背面から	3–4
	STERI-CYCLE i160 LK / STERI-CYCLE i250 LK ループスルーガス供給:	3–5
	STERI-CYCLE i160 LK / i250 LK の安全装置	3–6
	装置内使用空間の環境	3–6
	温度:	3–6
	相対湿度:	3–6
	水質の推薦:	3–7
	プレフィルター	3–8
	HEPA フィルターとエアダクト	3–8
	CO2 供給:	3–9
	02 供給:	3–9
	N2 供給:	3-10
	ドアスイッチ	3-10
	センサー	3-11
	供給インターフェース	3-13
	標準インターフェース	3-13
	オプションインターフェース	3-14
	ガス接続端子:	3-15
	注意書き:	3-15
	USB インターフェース:	3-15
	インターフェース 4-20 mA:	3-15
	アラームコンタクト:	3-16
	電力接続:	3-16
	装置内使用空間のコンポーネント	3-17
	内側容器	3-17
	ガラス扉とオプションのガスシェード	3-17
	貯水槽	3-19
	インキュベーターへの給水	3-21
	ヒーターシステム	3-21
	裏側の装置開口部分	3-22
	棚システム	3-23
	電気機械ドアロック・キット	3-24
4	装置使用の開始	4-1
	装置を環境に順応させる	4-2
	装置内使用空間の処理をします	4–2
	注水マーク「MAX」及びプレフィルターの交換	4–3
	エアダクトの取り付け	4–4
	HEPA フィルター及び貯水槽カバーの取り付け	4–6
	棚システムを設置する	4–8
	プロフィールの取付け取外し	4–8
	取付金具の装着	4–9
	装置のラベリング	4–9
	分割インサートプレート(オプション)を STERI-CYCLE i250 LK	
	に取り付ける	4-10
	ガスを接続	4-10
	ガス耐圧ホースの取付け....................................	4-11
	ガス接続 ガスモニタリングなし	4-12
	CO2 接続 ガスモニタリングシステム (オプション)付き	4-13
	CO2 と O2/N2 のコンビ接続 ガスモニタリングシステム (オプション)付き	4-14

電力接続		
いい インダーン	⊥ ─ へ ど 按 枕 ・	
アラームコン	クトを接続:	
インターフェー	-ス 4-20mA に接続する∶	
運転		
装置の準備		
運転を開始		
インキュベー		
潘 転		
电 ノスイッナ	······································	
コントロール	ペイルと操作力法	
02/N2 調整な		
CO2/O2/N2 ¬	ルチガス(オプション):	
操作方法		
iCan™タッチス	クリーンの初期設定	
ループセンサ-	・のウォームアップ	
設定時のキー	ダン操作	
温度設定値を	定する	
(1) 設定値を	、2000	
002 改定値で	、た 7 ゆ · · · · · · · · · · · · · · · · · ·	
02 改た値で改	ニ 9	
日期人ダート	〖	
目動スター	·をオンにする	
目動スター	を停止する	
Steri-Run を呼	び出す	
ユーザーコン	'ィギュレーション	
設定 / セッ	トアップ	
イベント・I	グ	
オプション		
アイコン説明]	
キーロック	·オン / オフにする	
ソフトウェ	'のバージョン	
	- JL	
エラー表示		
エフ 衣示 …	べいトに対するリアクション	
ユリー 衣小・	ペントに対するワアフション	
「四四反休役」	:86 ノビノド	
「「一日夜の一世」		
エラーの原		
装置の利用停」	:	
装置の利用を	ら止する	
洗海レ消毒		
ルガC/月母・・		
元才	•••••••••••••••••••••••••••••••••••••••	
际栄ノロセム		
消毒と ster i-	un を準備する	
拭き消毒 / ス	[/] レー消毒	
Steri-Run 除鲜	Ξ -ド	
Steri-Run 🎼	染モードのプロセス:	
Steri-Run を	オンにする	
Steri-Run z	中断する	

Steri-Run のエラー中断 Steri-Run 終える Steri-run 除染モード中断後にドアを開ける	8–10 8–11 8–12
メンテナンス 検査とコントロール メンテナンスの間隔 温度キャリブレーションの準備 温度の整合を行う CO2 キャリブレーションを準備する. CO2 キャリブレーションを準備する. CO2 キャリブレーションを行う HEPA フィルター 交換. ガスダクトフィルター交換 装置安全装置の交換 ドアパッキンの交換.	. 9–1 . 9–1 . 9–2 . 9–2 . 9–3 . 9–5 . 9–6 . 9–7 . 9–9 9–10 9–10
廃棄処分	10–1 10–1
仕様 STERI-CYCLE i160 LK. STERI-CYCLE i160 LK . STERI-CYCLE i250 LK . STERI-CYCLE i250 LK .	11–1 11-2 11–4 11–6 11–8
添付物 データ通信 USB インターフェース ドライバ USB インターフェースをインストールする データコミュニケーションのコマンドシーケンス構造 記録の説明 一般パラメータの一覧(アドレス 0xxx) ー般パラメータを読み取る 一覧インキュベータ・パラメータ(アドレス n 2xxx) パラメータ(Basic)を読み取る パラメータ(basic)を読み取る パラメータ(内部機能)を読み取る パラメータ(内部機能)を読み取る パターン 記録構造 エラーメモリ: エラー表示の一覧は Hex コーデックにて エラー表示の一覧は Bit コーデックにて ー般装置ステータス、ループ温度、CO2: ループ 02 および水位: データログ構造 パターン 記録構造 データログ: イベント入力情報一覧パート I: イベント入力情報一覧パート I: コード例 データログ データログデータ確認の機能 データログデータ確認のコード例 プログラム STERI-CYCLE i160 LK / STERI-CYCLE i250 LK STERI-CYCLE i160 LK & STERI-CYCLE i250 LK	12–1 12–2 12–4 12–4 12–5 12–5 12–5 12–5 12–6 12–7 12–8 12–9 12–10 12–10 12–11 12–11 12–13 12–14 12–15 12–16 12–16 12–21 12–22 12–23
	Steri-Run のエラー中断. Steri-run 除染モード中断後にドアを開ける. メンテナンス 検査とコントロール. メンテナンスの間隔. 温度キャリブレーションの準備. 温度の整合を行う. C02 キャリブレーションを在き. C02 キャリブレーションを有う. HFPA フィルター交換. ガスダクトフィルター交換. 携案処分. 使用素材一覧: 仕様 STERI-CYCLE 160 LK. STERI-CYCLE 160 LK. STERI-CYCLE 160 LK. STERI-CYCLE 160 LK. STERI-CYCLE 1230 LK. STERI-CYCLE 1200 LK. STERI-CYCLE 1230 LK. STERI-CYCLE 1300 LK STERI-CYCLE 130 LK STERI-CYCLE 130 LK

13	装置ブック	13–1
14	Thermo Scientific のコンタクトデータ	14–1
14	Cell locker 用 6 ドア気密スクリーン 棚板の設置	14–1 14–2

×

义	2-1。装置の寸法	2–2
义	2-2。持上げ	2–3
义	2-3。装置を重ねる	2–4
义	2-4。枠とエレメント付の台座	2–5
义	2-5。アダプタープレートと下側の装置を固定	2–5
义	2-6。キャリアフレーム取り外し用ねじ	2–6
义	2-7。キャリアフレーム取り付け	2–6
义	2-8。足なしの可動式台座(i160 LK 運転時)	2-8
义	2-9。スタンドの強化	2–9
义	2-10。装置のサポートスタンドをカウンターウェイト(おもり)で固定	2–9
义	2-11。装置のサポートスタンドを後方の壁に固定	. 2–10
义	2-12。前面の装置スタンドを固定	. 2–10
义	3−1。STERI-CYCLE i160 LK / i250 LK 前方から	3–2
义	3-2。STERI-CYCLE i160 LK / i250 LK 背面から	3–4
义	3-3。STERI-CYCLE i160 LK / STERI-CYCLE i250 LK ループスルーガス供給	3–5
义	3-4。HEPA フィルターとエアボックス	3–8
义	3-5。エアダクト	3–9
义	3-6。ドアスイッチ	. 3–10
义	3-7。温度、02、C02、空気湿度用センサー	. 3–11
义	3-8。センサー 水位用	. 3–12
义	3-9。装置インターフェース (スイッチボックスの右側)	. 3–13
义	3-10。オプションの装置インターフェース (スイッチボックスの左側)	. 3–14
义	3-11。STERI-CYCLE i160 LK ガラス扉付き	. 3–18
义	3-12。STERI-CYCLE i160 LK 3 分割ガスシェード付き	. 3–19
义	3-13。貯水槽	. 3–20
义	3-14。給水・排水 貯水槽の排水バルブ	. 3–20
义	3-15。インキュベーターの排水・給水バルブ	. 3–21
义	3-16。給水用漏斗	. 3–21
义	3-17。裏側の装置開口部分	. 3–22
义	3-18。棚システムの部品	. 3–23
义	3-19。インキュベーター下部のドアロックレバーと緊急用ロック解除	. 3–24
义	4-1。MAX と書いてある注水マーク	4–3
义	4-2。注水マーク「MAX」及びプレフィルター	4–3
义	4-3。エアダクト 組立てる	4–4
义	4-4。エアダクト構成部品 STERI-CYCLE i160 LK	4-5

义	4-5。HEPA フィルター とエアボックス 組立てる	4-6
义	4-6。エアボックスを貯水槽カバーに乗せる	. 4–7
义	4-7。エアボックスを取付ける	. 4–7
义	4-8。取付け/取外し 棚システム	. 4–8
义	4-9。取付金具の装着	4–9
义	4-10。取付け 分割インサートプレート	4–10
义	4-11。ガス耐圧ホースの取付け	4–11
义	4-12。ガス接続 ガスモニタリングなし	4–12
义	4-13。CO2 接続 ガスモニタリングシステム (オプション)付き	4–13
义	4-14。CO2 と O2/N2 のコンビ接続 ガスモニタリングシステム (オプション)付き	4–14
义	4-15。電力接続	4–15
义	4-16。例 接続 アラームコンタクト	4–18
义	4-17。インターフェース 4-20 mA の接続レイアウト	4–19
义	5-1。貯水槽	. 5–2
义	5-2。インキュベーターの排水・給水バルブ	5–3
义	5-3。給水用漏斗	. 5–3
义	5-4。MAX と書いてある注水マーク	5-4
义	6-1。電力スイッチ	6-2
义	6-2。メインディスプレイ: タッチディスプレイ	. 6-3
义	6-3。iCan™タッチスクリーン ガス供給 02/N2 なし	. 6–4
义	6-4。iCan™ タッチスクリーン、コンビガス接続付き	. 6–5
义	6-5。メニュー構造一覧	6-6
义	6-6。ウォームアップの表示	. 6-7
义	6-7。温度表示と温度選択メニュー	. 6-8
义	6-8。温度設定値を設定する	6-9
义	6-9。CO2 表示と CO2 メニュー	. 6-9
义	6-10。CO2 設定値を設定する	6–10
义	6-11。02 表示箇所と 02 メニュー	6–11
义	6-12。02 設定値を設定する	6–11
义	6-13。自動スタートをオンにする	6–14
义	6-14。自動スタートのステータス表示	6–15
义	6-15。自動スタートを停止する	6–16
义	6-16。自動スタート停止後のエラー表示	6–16
义	6-17。ユーザーコンフィギュレーションメニュー	6–17
义	6-18。メニュー 設定 / セットアップ	6–18
义	6-19。キーロックのコードを変更する	6–18
义	6-20。キーロックのコードを変更する	6–19
义	6-21。選択メニュー 日付 / 時刻	6–20
义	6-22。日付を設定する	6–20
义	6-23。時刻を設定する	6–21

义	6-24。	ディスプレイの明るさを設定する	6-22
义	6-25。	キー音の音量を設定する	6-23
义	6-26。	選択メニュー インターフェース USB	6-24
义	6-27。	USB インターフェースのボーレートを設定する	6-24
义	6-28。	言語を設定する	6-25
义	6-29。	リマインダー機能を選択します。	6-26
义	6-30 _°	steri-run のインターバルを設定する	6-27
义	6-31。	選択メニュー データ記録	6-28
义	6-32°	イベントを表示する	6-28
义	6-33°	記録インターバルを設定する	6-29
义	6-34。	エラー表を見る	6-31
义	6-35。	選択メニュー オプション	6-32
义	6-36。	選択メニュー アラーム	6-32
义	6-37。	アラームリレーを設定する	6-33
义	6-38°	低湿度を設定する	6-34
义	6-39 _°	ガスシェードを設定する	6-35
义	6-40 _°	水位センサーを設定する	6-36
义	6-41。	アラームリレーを設定する	6-37
义	6-42。	02 調整をオン / オフにする	6-38
义	6-43°	HEPA- コンフィギュレーション	6-39
义	6-44°	HEPA フィルターをオン / オフにする	6-39
义	$6-45_{\circ}$	アイコン説明	6-40
义	$6-46_{\circ}$	ガスモニタリングのアイコン	6-41
义	647_{\circ}	キーロックをオン / オフにする	6-43
义	648_{\circ}	ソフトウェアのバージョン	6-43
义	649_{\circ}	CO2 濃度の経過表示 を呼び出す	6-44
义	650_{\circ}	CO2 濃度の経過表示 を呼び出す	6-44
义	6-51。	イベントエラー表示	6-46
义	6-52°	エラー表示 過温度	6-46
义	6-53°	エラー表示 過温度	6-47
义	7-1。#	給水・排水 貯水槽の排水バルブ	. 7–1
义	8-1. #	給水・排水 貯水槽の排水バルブ	. 8–2
义	8-2° f	HEPA フィルターとエアボックス	. 8–4
义	8-3° ;	エアダクト	. 8–5
义	8-4。 [除染モードのプロセス	. 8–8
义	8-5°	メニュー Steri-Run 操作方法	. 8–8
义	8-6. 3	Steri-Run をオンにする	. 8–9
义	8-7° S	Steri-Run を停止、中断する	8-10
义	8-8. 3	Steri-Run を終了	8-11
义	8-9 _° ·	インキュベーター下部のドアロックレバーと緊急用ロック解除	8-12

义	9-1 _°	温度キャリブレーションの準備	9–3
义	9-2°	温度表示と温度選択メニュー	9-4
义	9-3°	温度の整合を行う	9-4
义	$9-4_{\circ}$	ガスシェードの計測口	9-6
义	$9-5_{\circ}$	CO2 キャリブレーションを行う	9–7
义	9-6.	エアボックスを取り外す	9-8
义	9-7 _°	HEPA フィルターの取付け	9-9
义	9-8 _°	ガスダクトフィルター 取付け	9–10

前書き

一般事項

装置の情報及び記録

ID データ

装置 ID

装置名称: 型名:	CO ₂ インキュベータ Steri-cycLe i160 LK Steri-cycLe i250 LK
製品記録の分類 CD 媒体	50143971

品質監査の証明:

コンプライアンス:	CE コンプライアンス証明
検査マーク:	TÜV GS, cCSAus, EAC

利用者への指示

本書は、CO₂インキュベータ STERI-CYCLE i160 LK / STERI-CYCLE i250 LK の使用説明書です。本 CO₂インキュベータは、現在最新の技術をもって製造し配送前に機能に欠陥がないことを 検査しております。しかしそれでも、本装置は危険が伴う事があります。特に、使用者 に知識が不足している場合、また、指定された利用以外にご使用になった場合などです。 事故防止のためには、それゆえ、以下にお気を付けください:

- 本 CO₂ インキュベータは、然るべき教育を受けそして許可を得た者のみ、使用することができます。
- この装置使って作業する者のために、本使用説明書、有効な安全データシート、社内の衛生規約、ならびにオペレーターの該当する技術規則を基に、書面にて手順指示書を作成してください。特に:
 - 本装置に対し使用する除染処置および補助手段、
 - ・ガスや加圧ガス容器の使用の際に守るべき安全処置、
 - 事故発生時の対処法
- · 装置のメンテナンスは、充分な知識と許可のある専門職員のみ行うことができます。

説明書の有効性

翻訳用の原本

- ・本使用説明書の内容は、予告なく変更される場合があります。
- ・本取扱説明書の英語版と翻訳版とのあいだに不一致がある場合、英語版が効力を有 するものとします。
- ・本使用説明書は、装置の近くに保管し、使用の際に安全に関する注意やその他の重要な情報をすぐに確認ができる状態にしてください。

本仕様書の記載が充分でないと感じお問合せがあります場合は、ご自身の安全のため Thermo Scientific までご連絡ください。

保証

保証条件

Thermo Scientific は、 CO_2 インキュベータの安全と機能に関し、以下の条件以外でのご利用の場合は一切責任を負いかねますので、予めご了承ください(2年間):

- ・本装置は、規定された利用目的にのみ使用され、本使用説明書に記載された方法で 利用ならびに維持すること、
- ・装置を改造しないこと、
- Thermo Scientific において認められている純正のスペアパーツおよび付属品を使用 すること、
- ・ 検査及びメンテナンスは、定められて機関で実施すること。

Cell locker およびすべての付属品には1年間の保証がついています。保証期間は、装置 が発注者に引き渡された時点から開始となります。

安全上のご注意とマークの説明

本使用説明書における安全上のご注意とマークの説明



留意 ご利用上のヒントや便利な情報を示しています。

安全上の注意を示すマーク:



保護手袋を着用!



保護めがねを着用!



危険な液体!



感電注意!



表面熱い!



発火の危険!



爆発の危険!



窒息の危険!

装置に記載のあるマーク

CE コンプライアンス証明: EU の要綱に沿ったコンプライアンスの証明です



TÜV によって検査された安全性



アメリカ合衆国・カナダの検査マーク



EAC 適合証明は、ユーラシア関税同盟(ロシア、カザフスタン、白ロシア)の全技術的 規定に適合することを証明するものです。



使用説明書に注意!



表面熱い!

装置の利用目的

規定の利用目的

本 CO₂ インキュベーターは、細胞の処理と培養用に考えられておりますが、医療及び体 外診断用ではありません。装置の利用場所は、本使用目的のため、以下のパラメータに よって的確に制御し、管理された生理学的環境条件を整えること:

- ・ 温度
- · CO₂含有量
- · 0₂/N₂含有量
- · 相対湿度

本 CO₂ インキュベーターの主な設置利用分野は、以下の通りです。

· 安全レベルL1、L2、L3の実験を行う細胞学や生物工学のラボ

- DIN EN 12128 に従った医学や微生物学のラボ
- ・クリニックや病院の研究ラボ

本装置は、専門的な使用にのみ適しており、教育を受けた職員のみが運転使用を許されています。

利用目的に反した使用

安全レベルL1、L2、L3の規定に沿わない細胞と組織培養を行ってはいけません。また、 以下の組織、素材、液体をサンプルとして使用してはいけません:

- ・ 容易に発火および爆発する危険性のある物、
- ・発する蒸気が空気に触れることで発火や爆発するような混合ガスとなる物、
- ・毒が発生する物。
- Cell locker は医療機器として使用しないでください。

規定と要綱

本装置は、以下の規定と要綱の安全要求に沿っています:

- · 定電圧指令 2014/35/EG
- IEC 61010-1:2010+Korr.2011、電気測定、制御、調整及びラボ装置に対する安全要 求事項、第1部:一般要求事項
- IEC 61010-2-010:2003、電気測定、制御、調整及びラボ装置に対する安全要求事項、
 第2部-010:素材を熱するラボ装置に対する特別要求事項
- · EMV 要綱 2014/30/EU
- IEC 61326-1:2012、電気測定、制御、調整及びラボ装置に対する安全要求事項、第 1部:一般要求事項
- ・ FCC 47、第 15 部、 § 15.107、EMV 規制 FCC
- · FCC 47、第 15 部、 § 15.109、EMV 規制 FCC

外国においては、当地の国内規約に拘束されます。

US (FCC)

"NOTE: This equipment has been tested and found to comply with the limits for a Class A digital device, pursuant to part 15 of the FCC Rules. These limits are designed to provide reasonable protection against harmful interference when the equipment is operated in a commercial environment. This equipment generates, uses, and can radiate radio frequency energy and, if not installed and used in accordance with the instruction manual, may cause harmful interference to radio communications. Operation of this equipment in a residential area is likely to cause harmful interference in which case the user will be required to correct the interference at his own expense."

Canada (ICES-001)

"This ISM device complies with Canadian ICES-001. Cet appareil ISM est conforme à la norme NMB-001 du Canada."

ガスに対する安全上の注意

留意 取付け:

 CO_2 や O_2/N_2 が入っている供給管ならびに加圧ガス容器、ボトル、収容器の扱いは、 専門職員によって適切な道具を使用したうえで行ってください。

留意

CO₂ 及び O₂/N₂ の作業場限界値を超えないようご注意ください。 ドイツ連邦共和国においては、TRGS 900 を守ることをお勧めします。その他の国に関 しては、これとは異なる限界値が有効となっている可能性があります。その国内で各 国特有の作業場限界値を遵守してください。

職員への指示:

装置 CO₂ 供給の装置を使用する職員は、この仕事を担当する前に CO₂ の扱いについて教育 を受けること:

- ・加圧ガス容器およびガス供給の適切な使用方法、
- · CO₂供給管における損失や血管の報告義務、
- · 事故や故障の際に行うべき処置。

これらの教育は、適度な期間を開けて繰り返し行ってください。ガス供給会社の特別な 取扱指示においても、教育内容に加えること。

二酸化炭素における安全上の注意(CO₂)

窒息の危険!

CO₂は、健康を損なう可能性のあるガスとしてレベル付けされています。それゆえ、本 CO₂インキュベータの投入および使用中において、以下の安全対策を守ってください:



二酸化炭素(CO₂)が大量に空気中に放出されると、窒素の危険があります。 CO₂流出の際は、すぐに安全対策を実行してください!

- · 該当の部屋からすぐに退出し、出入り口の確保をしてください!
- ・ 警備員もしくは消防署に連絡してください!

酸素における安全上の注意(0₂)

02は、可燃性のガスであり、油分を含む素材と反応し爆発する可能性があります。

酸素爆発の危険!



酸素(02)は、油、油分、グリースと触れることで爆発の危険性があります。高濃度酸 素と油分を含む物質が接触する場合、爆発の危険が発生します! 操作および表示ディ スプレイフィルム!

・本装置のクリーニングには、油分を含まない洗浄剤のみご利用ください。

酸素設備の接続部や部品からは、油、油分、グリースが含まれている素材を離してお くこと!

発火の危険!



酸素(02)の流出は、発火の危険性が大きいのでご注意ください。酸素を使用する設備の周りでは、裸火を使用しないでください!

·酸素設備の近くでは喫煙しないでください。

酸素設備の部品は、熱の影響下にさらさないでください。

窒素における安全上の注意(N₂)

窒素は、空気と混ざりやすい物質です。高濃度の窒素は空気中の酸素量を減らしてしまいます。

窒息の危険!

窒素 (N_2) が大量に室内の空気に放出されると、酸素不足による窒息の危険が発生しま す。 N_2 流出時には、安全対策を実行してください!

· 該当の部屋からすぐに退出し、出入り口の確保をしてください!

· 警備員もしくは消防署に連絡してください!



コンテンツ

- ・「包装」上 頁 1-1
- ・「配送確認」上 頁 1-1
- ・「STERI-CYCLE i160 LK / i250 LK 通常装備の配送内容」上 頁 1-2
- 「STERI-CYCLE i160 LK / i250 LK 追加装備の配送内容」上 頁 1-2
- ・「STERI-CYCLE i250 LK 追加装備の配送内容」上 頁 1-3

包装

CO₂ インキュベータ STERI-CYCLE i160 LK / STERI-CYCLE i250 LK は、頑丈な梱包箱に入れてお届けいたします。梱包素材は全て、分別の上再利用が可能です:

- ・ 梱包箱 : 古紙
- · 合成物質発泡材:発泡スチロール(FCKWフリー)
- ・ 梱包シート: ポリエチレン
- ・ 梱包紐: ポリプロピレン
- ・ 基部: ポリプロピレン
- ・パレット:未加工の木材

配送確認

商品引き渡し後すぐに、以下を確認してください:

- ・配送物が全てあるか、
- ・装置の配送状態。

配送内容に不足がある場合や、運送時に装置への被害があった場合、特に湿気や水分に よる損害が確認できる場合は、配送業者及びのテクニカルサポートにすぐにご連絡くだ さい。

STERI-CYCLE i160 LK/ i250 LK 通常装備の配送内容

配送の装置部品	個数
貯水槽カバー	1
最高水位センサー	1
バッフル 背面	1
バッフルカバー	1
エアボックス、パッキン付	1
プレフィルター	1
インサートプレート	3
インサートプレート用プロフィール	4
インサートプレート用取付金具	6
カバーロゼット用の栓	1
電力ケーブル	1
接続ホースー式 CO ₂	1
使用説明書	1
排水用ホース付クイックリリースファスナー	1

STERI-CYCLE i160 LK/ i250 LK 追加装備の配送内容

酸素調整追加装備の配送内容	個数			
02 センサーヘッド	1			
0 ₂ 接続ホースー式	1			
ガスボトルモニター CO ₂ 追加装備の配送内容				
接続ホースー式 CO2 ガスボトルモニター	1			
ガスボトルモニター 0 ₂ /N ₂ 追加装備の配送内容				
接続ホースー式 0 ₂ /N ₂ ガスボトルモニター	1			
フィルター追加装備の配送内容				
HEPA フィルター または VOC フィルター	1			
インターフェース 420mA 追加装備の配送内容				
プラグ 420mA	1			

STERI-CYCLE i160 LK 追加装備の配送内容

コンポーネント	個数
3分割ガスシェード	1
6 分割ガスシェード	1
Cell Locker	6
給水用漏斗	1

STERI-CYCLE i250 LK 追加装備の配送内容

分割インサート付き 250 リッター装置追加装備の配送内容	個数
キャリアフレーム	3
分割インサートプレート	6
6 分割ガスシェード	1



コンテンツ

- 「周辺条件」上 頁 2-1
- ・「換気」上 頁 2-2
- ・「必要なスペース」上 頁 2-2
- ・「運送」上 頁 2-3
- ・「装置を重ねる」上 頁 2-3
- ·「追加装備、装備変更」上 頁 2-13

周辺条件

本装置は、記載された環境条件の整った設置場所でのみ、ご利用可能です:

要求事項:

- ・ 隙間風のなく乾燥した場所。
- ・ 隣の物は、必要最低限の間隔を必ず開けてください。参照 「必要なスペース」上 頁 2-2。
- ・装置使用部屋には、適切な換気扇がついていること。
- ・非可燃性の堅い面であること。
- ・装置の重さや装置の追加(特に統治を積み重ねる際)に対して耐久性があり振動のない土台(台座、ラボデスク)。
- · 本装置は、最高海抜 2000 m までの場所でご利用いただくことができます。
- インキュベーション温度が通常 37° C に維持されるよう、室内温度を
 +18° C ~ +34° C に設定してください。
- ・ 相対湿度は、最高 80 % まで。
- ・ 直射日光を避けてください。
- STERI-CYCLE i160 LK / STERI-CYCLE i250 LK の近くには、高熱を発生させるような装置を設置しないでください。

換気

CO₂/O₂/N₂の利用時は、圧力調整口から圧力が排出され、インキュベータ利用空間において若干の超過圧力が発生します。

圧力調整および運転中のガラス扉やガスシェードの開閉により、微量の CO₂/₀2/_N2 が室内 に放出されます。換気の際は、流出するガスを安全に屋外に排出するよう気を付ける必 要があります。

また、装置の継続運転の際、装置から発散されるエネルギーにより、室内の環境が変化 することがあります。

- STERI-CYCLE i160 LK / STERI-CYCLE i250 LK はそれゆえ、換気が充分な部屋に設置板ダウ必要があります。
- · 本装置は、換気のできない壁のくぼみ部分などには設置しないでください。
- ・ ラボに対する当地の要綱条件に合った技術換気設備、または、それに見合った性能 の高い換気設備をご利用ください。

必要なスペース



注意 緊急時オフ!

電力供給用のコンセント差込口は、緊急時の電力オフのため常にすぐにアクセスが可能な状態にしておいてください。電源接続は、使用者からはっきりとわかるところに、そして、すぐに手の届くところにすること。



図 2-1。 装置の寸法

留意 装置へのアクセス: ケアやメンテナンスのためにも、両脇および背面に充分なスペースを保ち、装置へは すぐにアクセスできるようお勧めします。回転軸(ヒンジ)開きのCell lockerと6 ドア気密スクリーンを備えたユニットは、ヒンジ側の壁から、31cm(12.2インチ)以 上間隔をあけておく必要があります。

運送

運送の際に、本装置も扉や背面のスイッチボックスなどのユニット部分を掴んで持たな いようご注意ください。



図 2-2。 持上げ

留意 持上げ: 本装置を持ち上げる際は、図に印の付いた場所のみを持つようにしてください。



注意 重量物です! 持ち上げるときは気をつけてください! 過度伸展や椎間円盤の損傷など、過負荷による怪我を避けるために、絶対に本イン キュベーターを一人では持ち上げようとしないでください。 物の落下による怪我を避けるため、インキュベーターを持ち上げるときは、安全靴等 の保護装備を着用すること。 指や手の(特に扉を閉めるときに挟んでしまっての)怪我や、インキュベーターの損 傷を避けるために、上記図に記されている持ち上げ点のみを使用してください。

装置を重ねる

STERI-CYCLE i160 LK / STERI-CYCLE i250 LK は、同じタイプの装置であれば最大 2 台まで重ね置き することが可能です。その際には、オプションの重ね置きアダプター(1/図 2-3)を間 に入れます。

オプションとして、装置を動かすための可動式台座(2/図 2-3)の配送が可能です。

代わりに、キャスターなしのその他のフレーム(4/図 2-3)を使用し、2 つのデバイス を積み重ねることもできます。

留意

積重ねアダプタプレートの取付け時、そして装置を重ねる際は、積重ねアダプタに同 封の取付け説明書をご覧ください。

留意 積重ねた装置の取扱い: キャスター付きの台座に乗せた装置は、床がまっすぐで段差がない限りは、同室内で あれば移動させて使用することができます。



図 2-3。 装置を重ねる

1。下に置く方の足付きの装置(3/図 2-3)を、積重ね器具(1/図 2-4)に乗せて、台座(2/図 2-4)上面に置いてください。



- 図 2-4。 枠とエレメント付の台座
- 2。スタッキング アダプター (7/図 2-5)の下面を、下側にくる装置の (8/図 2-5) 上面につけます。
- 3。スタッキング アダプター (9/ 図 2-5) の接続箇所 (7/ 図 2-5) の両側にある穴を、 下側の装置の背面にある穴に合わせて (8/ 図 2-5) 配置します。







- 図 2-5。 アダプタープレートと下側の装置を固定
- 4。2本のネジを使用して、接続タブ(9/図 2-5)付きスタッキングアダプター (7/図 2-5)を下部ユニット(8/図 2-5)の後方の壁に固定します。その後方の壁に 対する反対側の角に、同じようにして固定します。
- 5。上に重ねる装置の足を(6/図 2-5)アダプタープレートの上面にある(1/図 2-4)重 ね置きエレメントに置きます。
- 6。2本のネジを使用して、接続タブ(9/図 2-5)付きスタッキングアダプター (7/図 2-5)をアッパーユニット(10/図 2-5)の後方の壁に固定します。その後方の 壁に対する反対側の角に、同じようにして固定します。

サポートブラケットの取り付け



注意

付属品の設置ガイドに特に記載がない限り、すべての積み重ねたユニットにはサポー トブラケットを使用することをお勧めします。

1。上側の装置の背面上部にある4つのねじの取り外しは、図 2-6の矢印部分になりま す。



- 図 2-6。 キャリアフレーム取り外し用ねじ
- 2。 2 つの矩形継ぎ目板を固定するために、キャリアフレームの両端(4/図 2-7)を約 90 度の角度に折り曲げます。
- 3。 4 つのねじはキャリアフレームを装置背面に固定するために使用します。継ぎ目板は 装置前面下方部にあります。
- 4。かならず、上のデバイスのデバイス スタンドが、積み重ねるアダプター上に正しく 位置合わせされていることを確認してください。
- 5。キャリアフレームは、負荷に耐え得る建物の一部に木製ねじと合わせ釘のような適切 な器具で固定してください。



図 2-7。 キャリアフレーム取り付け



注意 積み重ねた装置の横転および落下の危険性 装置の横転のおそれを防ぐためにも、キャリアフレームは必ず積み重ねた装置の負荷に 耐え得る壁に取り付けてください。設置は認定を受けた職員によってのみ行われます。 キャリアフレームを建物の一部に固定する際は、負荷容量が少なくとも 25kg ある適 切なねじおよび合わせ釘を使用してください。



注意 重ねた装置の運送! 重ね置きエレメントは、接続具ではありません。キャスター付きの台座に乗せた装置 は、床がまっすぐで段差がない限りは、同室内であれば移動させて使用することがで きます。

留意 可動式台座の固定: 装置が可動式台座に設置されている場合、インキュベータが運転されている際にキャ スターのストッパー(使用する場合)がかかっていること、また、安全のためキャス ターが前向きになっているようにしてください。

重ねた装置の運転におけるコンデンセートの形成: 基本的に、*STERI-CYCLE i160 LK* 及び *i250 LK* の型式を積重ねて運転する場合は、断熱のためにアダプタプレートを使用する必要があります。 積重ねた装置を 28 °C 以上の周辺温度において運転する際に、下側の装置が Ster i-run 除染中モード中の場合、上側の装置においてオーバーヒートのエラーが発 生します。その際、上側の装置においてコンデンセートが形成されることがありま す。

重ね置きヴァリエーション

重ね置き可能な組み合わせ		積み重ね下部			
		HERAcell Vios 160i LK	HERAcell Vios 250i LK	HERAcell 150i	HERAcell 240i
		Steri-Cycle i160 LK	Steri-Cycle i250 LK		
	HERAcell Vios 160i LK Stori Cvolo	50148171	50154522	50148172	
積み重ね上部	i160 LK				
	HERAcell Vios 250i LK		50154522		50148175
	Steri-Cycle i250 LK				

詳細は設置説明書の積重ねアダプタ項をご参照ください。

2 装置の設置 サポートブラケットの取り付け

可動式台座の取り付け





図 2-8。 足なしの可動式台座(i160 LK 運転時)



注意安定性の保証 装置作動中は台座の脚輪がロックされ、前方に向いていることを確認してください。

- · 台座は脚輪上部にあるレバー (3/図 2-8) でロックしてください。
- デバイス スタンドを使用してユニットを位置合わせします(1/図 2-8)は4つの重ね置きプレート(2/図 2-8)の上に正確に合わせてください。
 運送の際に、本装置も扉や背面のスイッチボックスなどのユニット部分を掴んで持たないようご注意ください。
- · 下部ユニットの下背面の端にある 4 個のスクリュー(5/図 2-8)を取り外します。
- ・ 背面のブラケット(4/図 2-8)を下部ユニットの背面の壁およびサポート スタンド にねじ止めします。

可動式台座と足(i160 LK と i250 LK での例)

納品内容



図 2-9。 スタンドの強化

サポート スタンド上にユニットを配置します。

i160 LK

- M6x20 スクリュー(5/図 2-10)2 個を使用し、サポート スタンド(1/図 2-10)上に追加のウェイト(2/図 2-10)をねじ止めします。
- 下部ユニット(6/図 2-10)の下背面の端にある 4 個のスクリュー(3/図 2-10)を 取り外します。
- ・ M6x20 スクリュー (5/ 図 2-10) 2 個を使用し、接続タブ (4/ 図 2-10) 上を追加の ウェイト (2/ 図 2-10) にねじ止めします。
- 前に取り外したスクリュー(3/図 2-10)を使用して、接続タブ(4/図 2-10)を下 部ユニット(6/図 2-10)の背面の端にねじ止めします。



図 2-10。装置のサポートスタンドをカウンターウェイト(おもり)で固定

i250 LK

- 下部ユニット(2/図 2-11)の下背面の端にある 4 個のスクリュー(1/図 2-11)を 取り外します。
- 前に取り外したスクリューを使用して、接続タブ(4/図 2-11)を背面の下部ユニット(2/図 2-11)にねじ止めします。提供されているスクリューを使用して、接続タブ(4/図 2-11)をサポートスタンド(3/図 2-11)にねじ止めします。



- 図 2-11。装置のサポートスタンドを後方の壁に固定
- 前面ブラケット(1/図 2-12)をデバイス スタンド(4/図 2-12)の上部品
 (5/図 2-12)および下部品(6/図 2-12)の間に入れます。
- · 4 個のスクリューを使用して、前面ブラケット(1/図 2-12)をサポート スタンド (2/図 2-12)にねじ止めします。



図 2-12。前面の装置スタンドを固定

- ・ 台座(2/図 2-12)に搭載された装置を望ましい設置場所へ注意深く移動させます。
- ・ 台座 (2/図 2-12) が水平に位置するまでレンチで足 (3/図 2-12) を留めつけます。

留意

台座 (2/図 2-12) 上の装置を稼動させる前に、足 (3/図 2-12) をレンチでしっかり と留めつけてください。



注意安定性の保証

装置がお望みの設置場所へ置いた後、レンチで足の(図 2-9)ねじが取り外された ことを確認してください。



留意 装置は図に示された持ち上げ位置からのみ持ち上げてください。



注意 重量物です! 持ち上げるときは気をつけてください! 過度伸展や椎間円盤の損傷など、過負荷による怪我を避けるために、絶対に本イン キュベーターを一人では持ち上げようとしないでください。 物の落下による怪我を避けるため、インキュベーターを持ち上げるときは、安全靴等 の保護装備を着用すること。指や手の(特に扉を閉めるときに挟んでしまっての)怪 我や、インキュベーターの損傷を避けるために、上記図に記されている持ち上げ点の みを使用してください。



注意 可動式台座上の装置の横転。



装置の稼働中は細心の注意が必要です。不注意な装置の減速または加速は可動式台座 からの横転をひき起こすおそれがあります。装置ドアを開けたまま稼動させないでく ださい。可動式台座は実験室内の設定場所近辺にのみ使用されるもので、運搬移動を 目的としたものではありません。

パートナ ンバー	サポート スタンド オプション	説明
50165306		HERAcell VIOS 160i LKと Steri-Cycle i160 LK の可動 式台座
50163093		HERAcell VIOS 160i LKと Steri-Cycle i160 LK の可 動式台座
50163094		HERAcell VIOS 250i LK と Steri-Cycle i250iLK の可 動式台座
50145435		HERAcell VIOS 160i LK および Steri-Cycle i160 LK 用ダブルチャンバー・サポートフレーム、高さ 200mm (キャスターなし)
50145436		HERAcell VIOS 250i LK および Steri-Cycle i250 LK 用シングルチャンバー・サポートフレーム、高さ 780mm(キャスターなし)
50149102		HERAcell VIOS 160i LK および Steri-Cycle i160 LK 用ダブルチャンバー・サポートフレーム、高さ 200mm (キャスターなし)
50149125		HERAcell VIOS 250i LK および Steri-Cycle i250 LK 用シングルチャンバー・サポートフレーム、高さ 780mm(キャスターなし)
追加装備、装備変更

通常版の装備追加は、以下の部品において行うことが可能です:

STERI-CYCLE i160 LK

- ・3分割ガスシェード(ガスシェードをガラス扉の代わりに取付け)
- ・外扉とガラス扉のストッパー変更、
- · Cell locker と6ドア気密スクリーン、
- 施錠可能な外扉、
- ·赤外線(IR) CO₂ センサー(スタンダード搭載の WLD センサーに代えて)、
- · 計測データインターフェース 4-20 mA
- ・ガスボトルモニター

STERI-CYCLE i250 LK

- ・3分割ガスシェード(ガスシェードをガラス扉の代わりに取付け)
- ・外扉とガラス扉のストッパー変更、
- 施錠可能な外扉、
- ・分割インサートプレート、
- ·赤外線(IR) CO₂ センサー(スタンダード搭載の WLD センサーに代えて)、
- · 計測データインターフェース 4-20 mA
- ・ガスボトルモニター

留意 装備変更:

追加装備ならびに装備変更は、Thermo Electron LED GmbH のテクニカルサービスのみ 行うことが可能です。 <mark>2 装置の設置</mark> 追加装備、装備変更

3



コンテンツ

- ・「STERI-CYCLE i160 LK / i250 LK 前方から」上 頁 3-2
- ・「STERI-CYCLE i160 LK / i250 LK 背面から」上 頁 3-4
- ・「STERI-CYCLE i160 LK / i250 LK の安全装置」上 頁 3-6
- ・「装置内使用空間の環境」上 頁 3-6
- ・「ドアスイッチ」上頁 3-10
- ・「センサー」上頁 3-11
- ・「供給インターフェース」上 頁 3-13
- ・「装置内使用空間のコンポーネント」上 頁 3-17

STERI-CYCLE i160 LK / i250 LK 前方から



図 3-1。 STERI-CYCLE i160 LK / i250 LK 前方から

- 1。圧力調整口 / 栓付きカバーロゼット
- 2。ガラス扉用ドアスイッチ
- 3。温度センサー
- 4。0₂ センサー(オプション)
- 5。ドアノブ ガラス扉またはガスシェード
- 6。ガラス扉またはガスシェード(オプション)
- 7。外扉

- 8。マグネット扉パッキン、交換可能
- 9。プレフィルター
- 10。排水口
- 11。足、高さ調整可能
- 12。パッキン、ガラス扉、交換可能
- 13。パッキンと HEPA フィルター付エアボックス
- 14. –
- 15。プロフィール
- 16。取付金具付きインサートプレート
- 17_°-
- 18。CO₂ センサー (IR)
- 19。ドアノブ 外扉
- 20。iCan[™] タッチスクリーン(コントロールパネル)
- 21。電力スイッチ
- 22。a = 電気機械式ドアロック / b ドア外側のロックフック(選択式ドアロック搭載の 装置のみ)

STERI-CYCLE i160 LK / i250 LK 背面から







図 3-2。 STERI-CYCLE i160 LK / i250 LK 背面から

- スイッチボックス、供給インターフェース付き、ガスコンビネーション接続用(オ プション)CO2 および O2/N2、ガスモニタリングなし
- 2。パターン: ガス接続 CO₂ および O₂/N₂、ガスモニタリングなし
- 3。0₂/N₂ ガスボトル
- 4。CO₂ ガスボトル
- 5。 第一 CO₂ ガス供給 A 、オプションのガスモニタリングの場合
- 6。 第二 CO₂ ガス供給 B 、オプションのガスモニタリングの場合
- 7。スイッチボックス、供給インターフェース付き、ガスコンビネーション接続用(オプション) CO_2 および O_2 /N₂、ガスモニタリング付き(オプション)
- 8。第一 0₂/N₂ ガス供給 A 、オプションのガスモニタリングの場合
- 9。 第二 0₂/N₂ ガス供給 B、オプションのガスモニタリングの場合
- 10。パターン: ガス接続 CO₂ および O₂/N₂、オプションのガスモニタリングの場合

11。分配接続、CO₂ ガス供給ループスル一用、装置3台との接続可能(装置タイプは無関係)

カバー:分配接続、0₂/N₂ガス供給ループスルー用、装置3台との接続可能、スイッチボックスの逆側

STERI-CYCLE i160 LK / STERI-CYCLE i250 LK ループスルーガス供給:





- スイッチボックス、供給インターフェース付き、ガスコンビネーション接続用(オ プション)CO2 および O2/N2 、ガスモニタリング付き(オプション)
- 2。分配接続、CO₂ ガス供給ループスルー用、装置3台との接続可能(装置タイプは無関係)
- 3。0₂/N₂ ガス供給のループスルー用分配接続 (カバー)、装置3台との接続可能 (装置 タイプは無関係)
- 4。耐圧ホース接続用 「字コネクタ
- 5。ガス供給ループスルー用耐圧ホース

- 6。パターン: ガス接続 CO₂ および O₂/N₂ ガスモニタリング付き (オプション)、CO₂ お よび O_{2 用}
- 7。*STERI-CYCLE i160 LK / STERI-CYCLE i250 LK* ループスルーガス供給; これには、ガスコンビ ネーション接続(オプション) CO_2 および O_2/N_2 が一つ必要です。

STERI-CYCLE i160 LK / i250 LK の安全装置

本装置は、以下の安全装置 が搭載されています:

- ・ドアスイッチは、ガラス扉の開扉時に、装置内使用空間ヒーターと CO₂/O₂/N₂のガス 流出を遮断します。
- オプションであるガスモニタリング使用時には、満タンのガスボトルへのガス供給
 ヘガス監視機が切り替わります。
- · 過温度保護機能機は、被害につながる過剰加熱の前におこる問題を防ぎます。
- ・圧力調整口は、装置内使用空間圧力の調整を行います。
- アラームリレーをオンにすると、聴覚視覚に訴える警告シグナルが、運転中のエ ラーを教えてくれます。

装置内使用空間の環境

インキュベータ装置内使用空間では、細胞と組織の処理と培養のため、特別な生物学的 環境条件を模擬実験されます。装置内使用空間の環境は、以下のファクターにより定め られます:

- ・温度
- 相対湿度
- ・ CO₂ 濃度
- · 0₂ 濃度(オプション)

温度:

装置が問題なく運転するために、の温度を最低でも18°Cに、インキュベーション温度 は最低でも室温より3°C高く設定して下さい。 ヒータィングシステムにより、インキュベーション温度を55°Cまでの間で調整しま す。独立ヒーター回路付きの内側容器ヒーターと別途の外扉ヒーターは、装置内使用空 間の壁と天井およびガラス扉やガスシェードにコンデンセートが形成されないようにし てくれます。

相対湿度:

装置内使用空間ヒーターは、水分の蒸発を促し、装置内使用空間内での一定の湿度を保 つ役割があります。装置運転には、推奨品質の濾過水の在庫を充分な量用意する必要が あります。

・ STERI-CYCLE i160 LK 及び STERI-CYCLE i250 LK の最大水量:30

水質の推薦:

装置が問題なく運転するために、貯水槽は、滅菌した蒸留水または適切に濾過した水で 満たしてください。水の伝導率は、1 ~ 20 μS/cm の範囲にあるようにしてください(電 子抵抗は 50 kOhmcm ~ 1 MOhmcm)。



注意保証の消滅!

塩素消毒された水道水や塩素を含んだ添加物を使用した場合は、保証が出来ません。 保証は、伝導率が1 ~ 20 μ S/cm の範囲外で電気抵抗が50 kOhmcm ~ 1 MOhmcm 外の純水(ウルトラピュアウォーター)を使用した場合にも保証出来ません。 ご質問は、Thermo Fisher Scientific テクニカルサービスまでお問い合わせください。



注意 加湿リザーバーへの水道水および超純水の使用禁止 加湿リザーバーに注入する水は、滅菌蒸留水、またはそれと同等のレベルで処理され た水を使用してください。使用する水の許容電気伝導率は、 1μ S/cm ~ 20μ S/cm です (抵抗率は、50K Ω ・cm ~ $1M\Omega$ ・cm でなければなりません)。pH は 7 ~ 9 でなければな りません。抵抗率が 18. $2M\Omega$ ・cm に等しいか、またはそれに近い 1 型超純水や脱イオ ン水はイオンをほとんど含まず、これを使用すると、内部の部品からイオンを溶解す るので、ステンレス、銅、ガラスに損傷を与えます。脱イオン水や 1 型水しか利用で きない場合は、弱アルカリ性の重炭酸ナトリウムの滅菌溶液を加えて pH を上げ、イ オンを加えます [84mg/l (1mmol/l) 推奨]。

注意 塩素系消毒剤の使用禁止

ステンレス鋼は腐食しにくいだけで、腐食していきます。多くの化学物質、特に塩素 や酸化活性を伴う誘導体は、ステンレス鋼に悪影響を及ぼします。

塩素系消毒剤や一定量の硫酸銅を消毒剤として含む水溶液の使用は、排水バルブのス テンレス製連結部や銅製ジョイントを損傷する可能性があるため、推奨されません。 内部を清掃するには、中性洗剤の水溶液でクリーニングし、水で洗い流して残留物を 除去することをお勧めします。内部の表面とパーツを希釈した第4級アンモニウム消 毒剤で拭きます。次に、70%アルコールで残っている微量の消毒液を拭き取ります。

通常の運転条件でインキュベーション温度が 37 °Cの場合、装置内使用空間相対湿度を約 93%に保ってください。

相対湿度が高いために培養容器にコンデンゼーションが形成されてしまう場合、装置内 使用空間の湿度を下げることができます。低湿度をオンにしている状態では、装置内使 用空間の相対湿度は約93%から約90%に下がります。変更をする場合は、調整に時間 がかかる可能性があります。培養容器のコンデンゼーションを効果的に避けるために、 継続的にこの設定にするようにしてください。

低湿度をオンにする点についての説明は、章「低湿度を設定する」上 頁 6-33 に記載が あります。

プレフィルター

貯水槽カバーの前側に、プレフィルター が付いています。プレフィルターは、シリコン 枠付きの二重金属メッシュでできており、オートクレーブ対応かつ耐熱です。Steri-Run 除染モード中は、プレフィルターは装置に付いたままにする必要があり、貯水槽への注 水の際に取り外します。

HEPA フィルターとエアダクト

貯水槽から装置内使用空間の方向への空気を流すことで、HEPA フィルターによって汚染 リスクを減少させることができます。フィルターは分離度 99.998 % で、粒子の大きさ 0.3 μm です(HEPA フィルター品質)。

HEPA フィルター (2/図 3-4) は、下からエアボックス (1/図 3-4) の中に取り付けます。 貯水槽カバー (2/図 4-6) の台の上にあるエアボックスを、エアダクトの方に引きます。



図 3-4。 HEPA フィルターとエアボックス

HEPA フィルター用アラームをオンにする方法は、章「HEPA フィルターをオン / オフにする:」上 頁 6-39 に別途記載があります。

エアダクトは、装置内使用空間内の背面(3/図 3-5)から上面に流れるファンからの空気 の流れを促し、温度が均等になる働きをします。エアダクトはそれと同時に、装置内使用 空間内に流れ込むプロセスガスを誘導し、ガスの混合が最適な状態となるようにします。



図 3-5。 エアダクト

エアダクトは、2つの金属プレートでできています:

- 1。上面チャンネル(1/図 3-5)
- 2。背面チャンネル(2/図 3-5)
- 3。装置内使用空間内の背面(3/図 3-5)

エアダクトと HEPA フィルターの取り外しには、工具は必要ありません。

CO₂供給:

細胞と組織が培養できる条件を確保するため、装置内使用空間内において CO₂を供給します。

重炭酸塩でバッファされた培養メディアの pH 値は、主に装置内使用空間内の CO₂ 含有量 で決まります。 装置内使用空間の CO₂ 含有量は、0 ~ 20 % の間で調整することができます。

供給される CO2の質は、以下の通りです:

· 純度 最低 99.5 %

・医療的ガス品質。

02供給:

CO₂ インキュベータを 21 % 以上の酸素で運転する場合、装置内使用空間に酸素を供給し ます (オプション)。 装置内使用空間の O₂ 含有量は、21 % ~ 90 % で調整できます。

高酸素濃度運転の場合は、前書きの章ガスに対する安全上の注意ページ7の発火防止の 注意をお読みください。 N₂ 供給:

運転中に酸素含有量を21%以下(空気酸素含有量)に下げる場合、装置内使用空間に窒素を供給します。装置内使用空間の02含有量は、センサーと関係なく調整できます。

ドアスイッチ



図 3-6。 ドアスイッチ

装置内使用空間開口部の上部エッヂに、ドアスイッチがあります。ガラス扉を開けるこ とでドアスイッチがオンになると、装置内使用空間へのガス供給とヒーティングが中断 します。コントロールパネルに通知メッセージが表示されます。 30 秒以上扉が開いた状態の場合、短いシグナル音が鳴ります。10 分以上扉が開いた状態 の場合、ブザー音が鳴り、アラームリレーがオンになります。 外扉は、ガラス扉を正しく閉めない限り、閉めることができません。

留意 ガスシェード: オプションのガスシェード付き装置の場合、上記記載のドアスイッチ機能は外扉を開 けた時点でオンとなります。

センサー

装置内使用空間背面には、ファンの羽根車とセンサーモジュールが搭載されています:

- ・装置内使用空間の温度および過温度保護機能を把握するセンサー(1/図 3-7)。
- · 0₂ センサー(オプション)は、装置内使用空間内の酸素含有量を把握するセンサー (2/図 3-7)。
- CO₂ センサーは、装置内使用空間内の CO₂ 含有量を把握するセンサー(3/図 3-7)。
 装備によって、この位置に WLD センサーまたは IR センサー(オプション)が取付けられます。
- rHセンサーは、装置内使用空間内の空気湿度を把握するセンサー(4/図 3-7)。rH
 センサー(オプション)は、WLDセンサーと湿度補填の役割を果たします。IRセンサーを使用しての運転の場合は、これはありません。



- 図 3-7。 温度、0₂、C0₂、空気湿度用センサー
 - ・水位センサー(1/図 3-8)は、貯水槽(2/図 3-8)に水を足す必要がある場合に警告をだします。貯水槽の水が0.50まで減ったら、ディスプレイのrH欄にエラー rH 水の「エラー表示」上頁 6-45 不足 という表示が出ます(も参照)。



図 3-8。 センサー 水位用

加えて、注ぎ足し時期がわかるよう、最大水位のメカニック / 視覚的表示もついています(参照 / 図 3-8)。

装置内使用空間温度センサー および CO₂ センサーとオプションの O₂ センサーは、装置コントロールシステムの一部です。その中で、計測値が設定値と比較されます。コントロールシステムはこのデータを基準とし、ヒーティングと CO₂/O₂/N₂ ガス供給を調整します。ファンにより、ガスと湿度を含んだ空気が混ざり、装置内使用空間内の温度が均等になります。

過温度保護機能は、工場側で既にプログラミングされ、訓練されたサービス員のみ変更 を行うことができます。過温度保護機能は、培養を加熱の超過から防ぎます。 設定温度から1°C以上超えた場合、過温度保護機能がオンになり、装置内使用空間温 度が自動的に設定値まで下げられます。すなわち、インキュベーション運転はエラー時 も継続します。過温度保護機能が起動する度に、同時に視覚警告シグナルを発します。過 温度保護機能がオンになると:

・エラー表示(温度)実測値が高すぎる)とブザー音が発動し、

アラームリレーがオンになります。

エラー表示が反応すると、過温度保護機能がオンになったという印として、ディスプレ イに過温度のアイコンが表示され、温度表示部分が赤くなります。

供給インターフェース

標準インターフェース



図 3-9。 装置インターフェース (スイッチボックスの右側)

供給接続は全て装置背面の供給インターフェース (スイッチボックス) に搭載されてい ます。

(図 3-9) スイッチボックス の右側に、基本装備およびオプション装備の幾つかがついています:

- 1。オプションのガスモニタリング運転および他装置への CO₂ ガス供給ループスルー用 の内部ガス分配接続端子 4 つ
- 2。0₂ 接続ノズル(オプションのガスモニタリング付き CO₂ および O₂/N₂ ガス供給の場合 は、なし)
- 3。CO₂ 接続ノズル
- 4。注意
- 5。USB インターフェース
- 6。アラームコンタクト
- 7。電力接続

オプションインターフェース



図 3-10。オプションの装置インターフェース (スイッチボックスの左側)

左側の接続端子は(図 3-10)該当のオプション搭載の装置にのみ搭載されています:

- 1。オプションのガスモニタリング運転および他装置への 0₂ ガス供給ループスルーの内 部ガス分配接続端子 4 つ
- 2。インターフェース 4...20 mA (オプション)
- 3。0₂ 接続ノズル(オプション:オプションのガスモニタリング付き _{C0}2 _{および 0}2/N2 ガ ス供給の場合のみ)

ガス接続端子:

装置とガス供給設備の間のガス供給は、同封の接続ホースにて行います。CO₂ および O₂/N₂ は、接続ノズル(2 と 3/ 図 3-9 および 3/ 図 3-10)。 にて装置供給されます。プロセスガスは全て、最低 0.8 ~ 最高 1.0bar の範囲内で供給 される必要があります。この値は、装置に既に設定され、変更が不可能な物です。 装置内使用空間へのガス流入前に、分離度 99.998 %、粒子の大きさ 0.3 μm(HEPA フィル ター品質)のガスダクトフィルターを通します。 ガスコンビネーション接続(オプション)、ガスモニタリングシステム(オプション) なしは、図のとおりです。

注意書き:

図(4/図 3-9)は、ガス供給、アラームコンタクト接続端子のレイアウト、装置の安全処置についての記載です。

USB インターフェース:

USB インターフェース (5/図 3-9)を使用し、インキュベータをパソコンを接続することができます。USB 1.1 / USB 2.0 / USB 3.0 フルスピードコンバーチブルの接続により、すぐに (一時的な利用の場合も)重要なパラメータ (温度、 $CO_2/O_2/N_2$ 濃度、エラーコード等)にアクセスすることができます。

インターフェース 4-20 mA:

オプションのインターフェース 4-20 mA (2/ 図 3-10) は、デジタルに表示されている温度、CO₂ 濃度 および O₂ 濃度(オプション)の実測値を、出力電力各々 4-20 mA に変換します。ここでは 4 回路 D/A 変換器に関連してマイクロプロセッサが搭載されています。 D/A 変換器は、回路につき 16 bit の解像度で、65536 ステップに相当します。

インターフェースは、使用者側で 4-20 mA 入力の様々な外付け計測装置、例えば Thermo Scientific スマート Vue 無線モニタリングシステムなど(配送物には含まれていません)と接続が可能です。

出カシグナルは、以下の表のとおりです:

回路番号	出カパラメータ (実測値)	解像度	計測範囲	出力 シグナル	
1	温度	0.08 MA/° C	0200° C	420mA	
2	C02 濃度	0.8 mA/%	020%C0 ₂	420mA	
3	0 ₂ 濃度 *	0.16 mA/%	0100%0 ₂	420mA	
4	フリー				
* がついているパラメータは、該当オプション搭載時のみ出力されます。					

機能の後に計測値:

$$MW = I_0 + I \times \frac{I_{max} - I_0}{MB}$$

I₀ = 4m A、 MW = 計測値(°C、%、rH%) I_{max} = 20m A、MB = 計測範囲

出カシグナルの意味は、以下の表のとおりです:

電力	意味。
4 mA	計測値は、計測範囲の最低値と同様もしくはそれより小さい値です。
20 mA	計測値は、計測範囲の最大値と同様もしくはそれより大きい値です。
2 mA	本装置は全機能作動する状態ですが、有効な値がありません(例え ば、オプションのセンサーが搭載されていない、本装置がまだ ウォームアップ段階であるなど)。
0 mA	エラー

4 つの 4-20 mA 出力シグナルは、差し込み可能な8 ピンの接続ブロック(3.5 mm)に通じ ます。該当の組み合わせコネクタは、配送内容に含まれています。ピンのレイアウトは、 以下の表のとおりです。

ピン番号	回路名称
1	回路 1: 温度
2	マス回路 1
3	回路 2:CO ₂
4	マス回路 2
5	回路 3:0 ₂
6	マス回路 3
7	回路 4: フリー
8	マス回路 4

アラームコンタクト:

本装置は、使用者側で外付けの通信システム(例えば電話、建物ガイダンスシステム、 視覚もしくは聴覚的アラーム機)に接続が可能です。

加えてポテンシャルフリーのアラームコンタクトが装置に搭載されています。このコン タクトは、装置背面のスイッチボックスに通じます図 3-9(6/)。

留意 アラームコンタクト: アラームコンタクトは、ループで発生したエラーの場合にオンになります(参照エ ラー表示の章)。

電力接続:

装置の電力接続は、ケーブル mit IEC プラグケーブルをスイッチボックス の IEC コンセントの差込口(7/図 3-9)に差し込むことで行われます。電源接続は、使用者からはっきりとわかるところに、そして、すぐに手の届くところにすること。

装置内使用空間のコンポーネント

インキュベーター装置内使用空間は、技術的対策により水分凝縮が起こらないよう、また、装置内使用空間に直接組み込まれた HEPA フィルターシステムに妨げとなるような汚染が発生しないように設計されています。装置内使用空間の HEPA フィルターシステムは、湿度を上げるために使用する水を保護し、クリーンルームに似た質の空気(クリーンルームクラス ISO 5)を作り出します。その際に培養菌の保管エリアに影響を及ぼすことは全くありません。

内側容器

インキュベータ装置内使用空間は、インキュベーション運転を妨げる汚染が起こらない よう、設計されています。これは、コンデンセートの形成阻止および装置内使用空間に 組み込まれた HEPA フィルタリングシステムにより可能となります。このシステムによ り、菌の培養に使用できる面での切断なしで湿度調整に使用する貯水を保護し、そして ISO 5 に沿ったクリーンルームの質に空気を整えることができます。

内側容器の素材

標準バージョンは、ステンレス鋼または銅製の内側容器一つが搭載されています。

内側容器の素材によって、エアダクトや棚システムなどの装置内使用空間のコンポーネントの素材も、同様のステンレス鋼または銅素材となります。

HEPA フィルター用のエアボックスは耐熱の合成物質で作られており、Steri-Run 除染 モード作動の際にも使用する必要があります。



注意 HEPA フィルターのは、60°C までの耐熱で、オートクレーブではないため、 Ster i-Run 除染の前に外してください。

留意 銅部分の酸化:
 熱および湿度の影響で、内容容器の銅素材は酸化します。銅部分それゆえ、装置検査のためのテスト運転の時から変色が始まります。
 酸化層は、定期的な洗浄の際には取り除くべきでありません。銅素材の抗菌効果は、これによるものだからです。

棚システムの棚システム、エアボックス、エアダクト、貯水槽の蓋は、道具を使用する ことなく簡単に取り外すことができますので、装置の洗浄と手動で行う消毒は、処理が され面積も少ない内側容器のみとなります。

ガラス扉とオプションのガスシェード

標準装備として、装置 **STERI-CYCLE i160 LK** および **STERI-CYCLE i250 LK** には安全効果ガラスの扉が一つ搭載されています。



図 3-11。STERI-CYCLE i160 LK ガラス扉付き

オプションのガスシェードが装備されている装置は、サンプルを取る開口部がかなり小 さくなっているために、汚染のリスクが4倍も低くなります。また、以下インキュベー ションパラメーターのインターバルも短く済みます。

- · 装置内使用空間温度、
- · CO₂濃度
- · 0₂/N₂ 濃度、
- · 相対湿度

ガスシェードは、両型式のオプション装備としてご注文が可能です:

- · STERI-CYCLE i160 LK: 3 分割ガスシェード (参照 図 3-12)、
- ・ *STERI-CYCLE i250 LK*: 6 分割ガスシェード
- ・オプションの Cell locker と 6 ドア気密スクリーン付録参照。



図 3-12。STERI-CYCLE i160 LK 3 分割ガスシェード付き

ガスシェード(2/図 3-12)のガラス扉には、専用のピンが付いています(1/図 3-12)。

貯水槽

貯水槽は、内側容器底に組み込まれており、装置内使用空間とはカバーで区切られてい ます。排水口(2/図 3-13)は、貯水槽の前側にあり、装置前側にある排水バルブ(3/ 図 3-13)を通して速やかな排水が行われます。

貯水槽カバーの前側には、プレフィルター(4/)図 3-13 が取り付けられています。プレフィルターは、シリコン枠付きの二重金属メッシュでできており、オートクレーブ対応かつ耐熱です。Steri-Run 除染モード中は、プレフィルターは装置に付いたままにする必要があり、貯水槽への注水の際に取り外します。Steri-Run 除染モード中は、プレフィルターは装置に付いたままにする必要があり、貯水槽への注水の際に取り外します。

6ドア気密スクリーンで給水する場合の補足説明付録参照。



図 3-13。貯水槽

貯水槽は、章「センサー」上頁 3-11 に記載のある水位センサーで監視されています。

注水がどこまで可能かを示す、MAX と書いてある注水マーク(1/図 3-13)が貯水槽にかかっています。貯水槽の最大注水量は3 Qです。

インキュベーション中に水の交換を行う場合に、装置内使用空間のエラー発生をできる だけ抑えるため、本装置には前方に排水バルブが搭載されています。前方にある排水バ ルブに同封の排水ホースを付ければ、排水を即座に行うことができます。



図 3-14。給水・排水 貯水槽の排水バルブ

インキュベーターへの給水



図 3-15。インキュベーターの排水・給水バルブ インキュベーターの給水バルブから給水用漏斗で水を注ぐことができます。



図 3-16。給水用漏斗

上図のようにガラスドアのロックに給水用装置のフックを引っ掛けます。 ホースを排水・給水バルブに接続します。

連通管の原理に従って、給水用漏斗によってインキュベーターの水位の高低差が決定されます。給水用漏斗の水位線は、最小が 0.5L で、最大が 3.5L です。漏斗の最大容量は約 0.4L です。

- · インキュベーターの排水・給水バルブに漏斗を挿入します。
- 漏斗に少量の滅菌蒸留水を注いで、水位が安定するまで待ちます。これを数回繰り 返します。
- ・ 最大水位線に達するまで給水を繰り返します。

ヒーターシステム

装置内使用空間のヒーティングには、フラッとヒーティングシステムを使用します。発 熱体は、貯水槽上部のコンデンセート形成をできるだけ抑えるように位置しています。 装置の外扉と扉開口の範囲も同様にヒーティングされています。内部のガラス扉 / ガス シェードに熱を充てることで、コンデンセートの沈澱を防ぐことができます。 装置内使用空間内は、湿度が高くても常に見通せるようになっています。

裏側の装置開口部分



図 3-17。裏側の装置開口部分

栓で閉めることができるカバーロゼット(図 3-17042 mm)(1/図 3-17)を使うと、装置 内使用空間の線、ホース、追加センサーを移動させることができます。

圧力調整口(3/図 3-17)はカバーロゼット下、装置背面にあり、装置内空間と装置使用部屋の圧力の調整を行います。

留意

シリコンの磨耗を防ぐため、高温除染を始める前に、内側容器から必ずシリコンストッパーを取り外し、通路開口部の外側に差し込んでください。

留意 運転条件:

CO₂インキュベータ装置内使用空間にて補助装置を使用する場合、環境条件に関する 要求事項にご注意ください(表参照)。装置内使用空間にもたらされたエネルギーは、 温度の制御範囲開始に影響を及ぼします。装置内使用空間に追加で熱源を付ける場合 は、コンデンセート形成(例えばガラス扉に)が発生する可能性があります。

もたらされたエネルギー	温度制御範囲の開始	
	一般	例: RT* = 21 ° C
0 W	RT + 3 ° C	24 ° C
5 W	RT + 6. 5 °C	27.5 °C
10 W	RT + 9. 5 ° C	30.5 °C
15 W	RT + 1 3 ° C	34 ° C
20 W	RT + 16 °C	37 ° C
*RT = 室温		

棚システム



図 3-18。棚システムの部品

(3/図 3-18) 棚システムのプロフィールは、42 mm 間隔でついています。取付金具(2/ 図 3-18) は必要な培養容器の大きさによって取付けることができます。インサートプ レート(1/図 3-18) には傾き保護装置とエクステンションリミットが追加されていま す。*STERI-CYCLE i250 LK*の装置の下側インサートプレートは、分割なしのものが付きます。 図 3-18 の左側をご覧ください。*STERI-CYCLE i160 LK*の装置には、充填口のある下側イン サートプレートが同封されています(4/図 3-18)。棚システムについては、章「装置使 用の開始」上 頁 4-1 に詳細が記載されています。

6ドア気密スクリーンの棚板を使用する場合の補足説明付録参照。

電気機械ドアロック・キット

電気機械ドアロック・キットは内蔵モーターで作動する回転式ロック(図3-19の詳細A)、内蔵の機械動作性緊急用始動レバー(詳細C)、そしてドア内側に取り付けられているロックフック(詳細B)からなります。

ドアロック・キットは Steri-run 除染モード中のインキュベーター装置内へのアクセス を防ぐ安全装置です。この安全装置はインキュベーター装置内温度が 65 °Cに達する、 あるいは超えると直ちにインキュベーターのドアをロックします。除染モードが終了し、 温度が 65 °C以下に低下すると、ドアロック・キットがドアロックを解除します。



図 3-19。インキュベーター下部のドアロックレバーと緊急用ロック解除

停電が起こった際、レバーは最後にあった位置に留まっています。電力供給の回復と除 染モードの再開後、温度が65°C以下に下がり次第、レバーがドアロックを自動的に解 除します。必要の際は、緊急用ロック解除レバーを側壁方向に引くことでいつでもドア ロックの解除ができます。緊急用ロック解除レバーは電力スイッチ横のインキュベー ター下部にあります。図 3-19の詳細Cは、レバーの位置を表示するために側壁から可視 化した図です。

装置使用の開始

コンテンツ

- ・「装置を環境に順応させる」上 頁 4-2
- ・「装置内使用空間の処理をします」上 頁 4-2
- ・「注水マーク「MAX」及びプレフィルターの交換」上 頁 4-3
- ・「エアダクトの取り付け」上 頁 4-4
- ・「棚システムを設置する」上 頁 4-8
- 「分割インサートプレート(オプション)を STERI-CYCLE i250 LK に取り付ける」上 頁 4-10
- ・「ガスを接続」上 頁 4-10
- ・「電力接続」上 頁 4-15
- ・「USB インターフェースを接続:」上 頁 4-16
- ・「アラームコンタクトを接続:」上 頁 4-17

4

装置を環境に順応させる

注意 装置を環境に順応させる!



装置使用の開始前には必ず本装置を環境に順応させる必要があります。

- ・本装置は、スイッチを入れる前に約2時間、予定装置使用部屋温度に設定されている装置使用部屋において設置しておく必要があります。
- ・装置ドアを開けます。

装置内使用空間の処理をします

CO₂ インキュベータは、殺菌状態で配送されません。それゆえ、本装置は運転前に除染す る必要があります。

装置内使用空間の以下の部品は、前もって洗浄し滅菌してください:

- ・プロフィール
- · 取付金具
- ・プレフィルター
- ・エアダクト
- ・エアボックス
- ・インサートプレート
- ・装置内使用空間の表面
- ガラス扉パッキン
- ・ ガラス扉 / ガスシェード

留意 除染:

装置における洗浄と消毒の詳細については、別途章にまとめてあります(参照 「洗 浄と消毒」上 頁 8-1)。



注意

スペア部品またはアクセサリー類とともに配送されているマニュアルの手順または技術データをよくお読みください。手順または技術データは、本マニュアルに記載の内容と異なることがあります。

注水マーク「MAX」及びプレフィルターの交換

注水マーク「MAX」及びプレフィルターは、工具なしで交換が可能です。

- 1。排水バルブ(3/図 4-2)のホースが取り外されていることを確認します。
- 2。 貯水槽前方にある排水口(2/図 4-2) が遮られていないかを確認します。これは、装置前面の排水バルブ(3/図 4-2) を通しての排水に必要となります。
- 3。注水マーク「MAX」(1/図 4-1)を貯水槽カバー(5/図 4-1)の溝にはめます。



図 4-1。 MAX と書いてある注水マーク

4。 プレフィルター (4/図 4-2) を貯水槽カバーに取り付けます。



図 4-2。 注水マーク「MAX」及びプレフィルター

エアダクトの取り付け

- エアダクトの上部(1/図 4-3)を、背部(2/図 4-3)を 図 4-3 に向けて、AからE のステップの通りに組立てます。その際に、図 4-3 のステップCにおいて、ポジ ション補助垂がダクト下部の該当の四角穴にかみ合うように気を付けてください。
- 2。背部の足のブリッジ(2/図 4-3)を、背面の両スタッドボルトに取付け、エアダクト を後ろに倒します。
- 3。上部側面のキーホール (ステップ G/ 図 4-3) を、装置内空間天井の留めピンとかみ 合うようにします。





図 4-3。 エアダクト 組立てる

図 4-4 は内側容器内に内蔵されたエアダクトを構成するパーツを示しています。



図 4-4。 エアダクト構成部品 STERI-CYCLE i160 LK

留意 STERI-CYCLE i160 LK の運転に際して、背面が固定されるまでエアダクトのパッキンを 取り外さないでください。背面のエアダクトは、ステンレス製の内側容器付の VIOS160-運転に必要とされる適切な気流に関する重要なパーツです。

HEPA フィルター及び貯水槽カバーの取り付け

HEPA フィルター (2/図 4-5) は、下からエアボックス (1/図 4-5) の中に取り付けます。 貯水槽カバー (2/図 4-6) の台の上にあるエアボックスを、エアダクトの方に引きます。



- 図 4-5。 HEPA フィルター とエアボックス 組立てる
- 1。HEPA フィルター(2/図 4-5)を平面上に置きます。
- 2。エアボックス(1/図 4-5)を左に傾け、左側の垂(3/図 4-5)と一緒に HEPA フィル ター(4/図 4-5)の該当の溝に差し込みます。
- 3。エアボックス右側の垂(5/図 4-5)を HEPA フィルターのかみ合い(6/図 4-5) に留 めます。
- 4。 パッキン (8/ 図 4-5) をエアボックスのパイプソケットの溝 (7/ 図 4-5) に取り付け、全体的にしっかりと押し込みます。

5。エアボックス(1/図 4-6)を貯水槽カバーの台(2/図 4-6)に乗せます。



- 図 4-6。 エアボックスを貯水槽カバーに乗せる
- 6。装置内使用空間底に貯水槽カバーを取り付けます。
- 7。貯水槽カバーの前側を上にあげ、背面方向にスライドさせます(1/図 4-7)。



- 図 4-7。 エアボックスを取付ける
- 8。貯水槽カバーを、止めがかかるまで背面方向に押します。カバーを槽の最終ポジショ ンまで、そしてエアボックスのパイプソケットをファン排気口までスライドさせま す。
- 9。貯水槽カバーの前方の縁を貯水槽に差し込んでいきます / 図 4-7)。その際、エア ボックスのノズルがファンの排気口に沿っていきます。

棚システムを設置する



図 4-8。 取付け / 取外し 棚システム

棚システムの設置には、道具は必要ありません。プロフィールはバネ圧で留めます。取 付金具をプロフィールにはめ、インサートプレートを取付金具にスライドします。

プロフィールの取付け取外し

プロフィールは、打ち出しにより横から入れて留めます。プロフィールの留バネは、その際上向きになるようにしておいてください。

- 1。下側の打ち出しのプロフィールはをつけ、装置内空間の壁にたたみ、双方の打ち出 しにプロフィールが乗っている状態にします。
- 2。上側の打ち出しの裏に、留バネをはさみます。
- 3。プロフィール取り外しには、垂部分の留バネを打ち出しから下側に引出し、プロ フィールを取り出します。

取付金具の装着



図 4-9。 取付金具の装着

- 1。取付金具をプロフィールのパーフォレーションにいれ、レイヤー棒が下向きになる ようにします。
- 2。取付金具の垂直部分双方が、プロフィールにあっているかを確認します。
- インサートプレートを入れる:
 - 1。傾き保護装置付きのインサートプレート(5/図 4-9)を、装置背面の方向を向けて 取付金具に押し入れます。傾き保護装置は、同時にインサートプレートのガイドと なる形です。
 - 2。 リミッター(6/図 4-9)が取付金具の上に来るよう、インサートプレートを軽く持ち 上げます。
 - 3。取付金具が双方の傾き保護装置問題なく入っているかを確認してください。

留意 STERI-CYCLE i160 LK の場合、充填装置付きのインサートプレートを、一番下の位置に入れてください。 6 ドア気密スクリーンの棚受けは長方形の穴にのみ設置してください。棚板の設置については付録を参照してください。

装置のラベリング

- 1。水準器を真ん中のインサートプレートもしくはロールレセプタクルの上に置きます。
- 2。調整可能な装置の足を同封のスパナー(幅 24 mm)で回し、インサートプレートが全 方向水平になるように調整します。装置の足の高さの町営は、左右前後可能です。

分割インサートプレート(オプション)を STERI-CYCLE i250 LK に取り付ける



図 4-10。取付け 分割インサートプレート

STERI-CYCLE i250 LK に 6 分割ガスシェードと 6 分割インサートプレートのオプションが付く 場合、標準装備品のホードの代わりに、横のサポートの取付金具にサポートフレームを 3 つ、各々にインサートプレート 2 枚ずつ取り付けてください。

- 両インサートプレート(1/図 4-10)をサポートフレーム(2/図 4-10)の上に取り 付けます(ステップA)。
- 2。サポートフレーム(2/図 4-10)を取付金具(3/図 4-10)にはめます(ステップB)。

ガスを接続

留意 ガス品質:

注意 過圧!

ガスは、以下の通りの質がなければいけません:

- ・ 純度 最低 99.5 %
- ・医療的ガス品質。



ガスは、装置に対し操作圧力最大で1 bar までにすること。ガスがこれより高圧で流入されると、装置内部のバルブがきちっと閉まらず、それゆえガス供給の調整に問題が出る危険があります。

ガス供給は、最低 0.8 bar 最高 1 bar に設定し、この圧が変更されないことを確実 にしてください!
ガス耐圧ホースの取付け



図 4-11。ガス耐圧ホースの取付け

ガス供給設備から装置へのガス供給は、同封のフレキシブルなガス耐圧ホースを使って 行います:

- 1。ガス耐圧ホースをガス供給設備の接続ノズルに差し込みます。
- 2。(3/図 4-11) ガスダクトフィルターの保護カバーを取り外します。
- 3。ホースクランプ(1/図 4-11)をガス耐圧ホース(2/図 4-11)にスライドさせ、ガス 耐圧ホース(4/図 4-11)をガスダクトフィルターの接続ノズル(5/図 4-11)に差し 込みます。
- 4。ガス耐圧ホースをホースクランプでガスダクトフィルター接続ノズルに固定します。

注意 圧力調整口

常に圧の調整が行われるよう、圧力調整口は排気システムに接続してはいけません。 圧力調整口のダクトの延長や迂回は行ってはいけません。

ガス接続 ガスモニタリングなし



図 4-12。ガス接続 ガスモニタリングなし

ガス供給設備から装置へのガス供給は、同封のフレキシブルなガス耐圧ホースを使って 行います:

CO₂ 接続:

 CO₂ 接続付きの装置の場合、ガス供給はガスダクトフィルター(2/図 4-12)に接続 パターン(1/図 4-12)のとおりに接続します。

 $CO_2 \ge O_2/N_2$ のコンビ接続 (オプション):

CO₂/O₂/N₂のコンビ接続の場合、供給管は接続パターン(3/図 4-12)のとおりに以下のように設置します:

· 0₂/N₂ 供給をガスダクトフィルター上部(4/図 4-12)、

· CO₂ 供給をガスダクトフィルター下部(5/図 4-12)。

留意 手動ガスアラーム:

ガスモニタリング (オプション)なしの装置の場合、自動ガスアラーム はありません。 そのため、ガスボトルのガス残量を毎日確認する必要があります。

CO₂ 接続 ガスモニタリングシステム (オプション) 付き

続パターン(7/図 4-13)通りに接続します。



図 4-13。CO₂ 接続 ガスモニタリングシステム (オプション)付き ガス供給設備から装置へのガス供給は、同封のフレキシブルなガス耐圧ホースを使って 行います。 オプションのガスモニタリングシステム搭載されている CO₂ 接続付きの装置の場合、接

CO₂ 接続:

プロセスガス CO₂ 似て運転する装置の場合、そして、オプションのガスモニタリングシステムが搭載されている場合、ガス供給は以下の通りに接続します:

- ・ ポジション 1/図 4-13:分配接続 CO₂ガス供給用、他の装置1台~3台に接続。ガス供給ループスルーには、それぞれの装置をガス耐圧ホースで繋いでください。
- ・ポジション 2/図 4-13:加圧ガス容器 B からのガス供給をガスモニタリングシステムの上部の接続端子につなげます。
- ・ポジション 3/図 4-13:加圧ガス容器 A からのガス供給をガスモニタリングシステムの下部の接続端子につなげます。

・ポジション4&5/図 4-13:ガスモニタリングシステムの排出は工場側で既に短いガス耐圧ホース(ポジション 6/図 4-13)を使ってガスダクトフィルターにつなげてあります。

留意 ガス供給のループスルー: ガス供給のループスルーでは、最大4の装置を CO_9 ガス供給1台で賄うことができます。

 $CO_2 \ge O_2/N_2$ のコンビ接続 ガスモニタリングシステム (オプション)付き



図 4-14。 $CO_2 \ge O_2/N_2$ のコンビ接続 ガスモニタリングシステム (オプション)付き

ガス供給設備から装置へのガス供給は、同封のフレキシブルなガス耐圧ホースを使って 行います。

CO₂/O₂/N₂ コンビ接続で、オプションのガスモニタリングシステムが搭載されている装置は、接続パターン(10/図 4-14)の通りに接続します。

0₂/N₂供給:

- ・ 0₂/N₂ ガス供給用分配接続、他の装置1台~3代との接続(カバー図 4-14)。ガス供給ループスルーには、それぞれの装置をガス耐圧ホースで繋いでください。
- ・ポジション 2/図 4-14:加圧ガス容器 Bからのガス供給をガスモニタリングシステムの接続端子につなげます。
- ・ポジション 3/図 4-14:加圧ガス容器 A からのガス供給をガスモニタリングシステムの接続端子につなげます。
- ・ポジション 1/図 4-14: ガスモニタリングシステムの排出は工場側で既に短いガス 耐圧ホースを使ってガスダクトフィルター(ポジション 8/図 4-14)につなげてあ ります。

CO₂供給:

- ・ポジション 4/図 4-14:分配接続 CO₂ ガス供給用 、他の装置1台~3台に接続。 ガス供給ループスルーには、それぞれの装置をガス耐圧ホースで繋いでください。
- ・ポジション 5/図 4-14:加圧ガス容器 B からのガス供給をガスモニタリングシステムの接続端子につなげます。
- ・ポジション 6/図 4-14:加圧ガス容器 A からのガス供給をガスモニタリングシステムの接続端子につなげます。
- ・ポジション 7/図 4-14: ガスモニタリングシステムの排出は工場側で既に短いガス
 耐圧ホースを使ってガスダクトフィルター につなげてあります。

留意 ガス供給のループスルー:

ガス供給のループスルーでは、最大4の装置(装置タイプ無関係)をそれぞれ CO_2 ガス供給もしくは O_2/N_2 ガス供給1台で賄うことができます。

電力接続



図 4-15。電力接続



注意 感電注意! 電気が通っている部分をふれると、感電し生命の危険となる場合があります。 プラグと電気ケーブルは、電力接続前に損傷がないかなどを確認してください。 損傷のある接続コンポーネントは、電力接続にしようしてはいけません!

本装置を、規定通りに設置しアース処理されている電力 にサーキットブレーカー B 16 を安全処置として使用し接続します。

電力接続:

- 1。電力に接続する前に、コンセントの差込口の電圧値が装置の電力スイッチ左側にあ るタイプシールの記載とあっているかを確認します。電圧(V)の記載と最大電気 (A)が合っていない場合、本装置は接続してはいけません。
- 2。IEC プラグを装置スイッチボックスのコンセントの差込口に差し込みます。
- 3。電力ケーブルのセーフティプラグを規定通りにアース処理し安全なコンセントの差込 ロに差し込みます。
- 4。電力ケーブルが引っ張られたり押し付けられたりしていないよう、確認してください。
- 5。電源コードの取り付け:電源接続は、使用者からはっきりとわかるところに、そして、すぐに手の届くところにすること。電源コードのコンセントプラグは、全電極の セパレータです。

USB インターフェースを接続:

装置にはシリーズとして USB インターフェースが搭載されています。パソコンへの接続 は、市場に出回っている USB 1 または USB 2 ケーブル(インキュベータ側: USB プラグ タイプ B、パソコン側: USB プラグ タイプ A)を使っておこないます。

USB インターフェースは、USB 1.1 に該当し、USB 2.0 と3.0 (フルスピード)とも適合 します。パソコンとインキュベータ間のデータのやり取りには、現行の Windows に既に 搭載されていない場合は、USB インターフェースを通しパソコンに該当のドライバをイン ストールします。ドライバのインストール方法は「添付物 データ通信」上 頁 12-1 に記 載してあります。

インターフェースの伝達速度は、必要であれば決められたボーレート(9,600、19,200、 38,400、57,600 ボー)の間で変更が可能です。ボーレートの設定方法は、「USB インター フェースのボーレートを設定する」上 頁 6-23 章に記載があります。

アラームコンタクトを接続:

留意 専門作業:

Thermo Scientificは、インストールが専門員によって行われた場合にのみ、本装置の安全性と機能を保証します。

装置の外部アラームシステムへの接続は、電気工学 / テレコミュニケーションの教育 を受け許可された専門員のみが行うことができます!

機能:

システムエラーおよび温度やガスループ内でのエラーの場合は、接続されている報告 / アラームシステムに警告が発信されます。ポテンシャルフリーコンタクト(1 チェンジャー)は、以下のサーキット用に形成されています:

アラームリレー:

サーキット	電圧	外部安全処置
サーキット電圧付	最大 250 V [~]	最大 6 A
SELV サーキット(参照 VDE 0100、 パート 410)	25 V ~	最大 2 A
	60 V =	最大 1 A
SELV E サーキット(参照 VDE 0100、 パート 410)	50 V ~	最大 1 A
	120 V =	最大 0.5 A
運転状況	コンタクト 3 - 2	コンタクト 3 - 1
運転状況 電力供給 アラーム 「オフ」	Х	0
運転状況 電力供給 アラーム 「オン」	0	X
エラー 電力供給ア ラーム 「オフ」	0	X
エラー 電力供給ア ラーム 「オン」	X	0
凡例:X: コンタクト	クローズ/0: コンタクト オ	ープン

留意 スイッチ構造: アラームリレーは、ループを通してのエラーの場合に常にオンになります(参照 「エラー表示」上 頁 6-45)。

接続例:



図 4-16。例 接続 アラームコンタクト

接続ケーブルに接続するプラグは、配送品には含まれていませんが、別途ご注文可能です。 通知システムの外部サーキットの操作圧力と安全処置の値は、以下の表のとおりです。

- 1。接続ケーブルの線それぞれは、電気接続プランの割り当て通りに挟みます。
- 2。外部通知システム用接続ケーブルのプラグを装置裏面のスイッチボックスにあるイン ターフェースに差し込みます。

インターフェース 4-20mA に接続する:

計測データインターフェース 4-20mA レイアウトは以下の図のとおりです:

2+1 0 28 GND 2 0 (07 + 3)0 CO2 GND 4 0 02/N2 + 50 02/N2 GND 6 0 7 0 8 0

4...20mA

図 4-17。インターフェース 4-20 mAの接続レイアウト

留意 專門作業:

Thermo Scientificは、インストールが専門員によって行われた場合にのみ、本装置の安全性と機能を保証します。

装置の外部アラームシステムへの接続は、電気工学 / テレコミュニケーションの教育 を受け許可された専門員のみが行うことができます!

計測データインターフェース 4-20 mA の装備追加は、Thermo Fisher Scientificのテクニカルサービスのみ行うことができます。

4 装置使用の開始 インターフェース 4-20mA に接続する:

5



コンテンツ

- ・「装置の準備」上 頁 5-1
- ・「運転を開始」上 頁 5-2

装置の準備

本装置は装置使用の開始における手順(参照 「装置使用の開始」上 頁 4-1)を行った場合にのみ、運転が許可されます。

装置チェック:

運転の前に以下の部品の装置チェックを確認してください

- ガスホースは接続フィルターにきちっとつないであり、ホースクランプで安全が保たれている。
- カバーロゼットがきちっと閉じている。
- ・ 圧力調整口には妨げるものがなく装置内使用空間に設置されている。
- ガラス扉のパッキンに損傷がない。
- ・ ガラス扉 / ガスシェードの計測口はきちっと閉じている。
- ・ 棚システムの部品が、安全に取付けられている。
- HEPA フィルター付きエアボックスとエアダクトが規定通りに設置されている。

装置内空間を除染する:



注意 Cell locker の使用耐熱温度! 「除染実行」の前に、Cell locker を取り出す必要があります。使用耐熱温度は最高 121°C (250°F) です。

 Steri-Run 除染モードを実行する(参照 「Steri-Run 除染モード」上 頁 8-6) また は、衛生要綱に沿って装置内使用空間を運転前に 除染します。 **留意**衛生要綱: 培養保護のために、装置内使用空間を社内で既定のある衛星要綱にそって運転の前に 洗浄し滅菌する必要があります。

水の貯蔵:参照「相対湿度:」上頁 3-6。

運転中に貯蔵水量の最低ラインを下回った場合は、水を注ぎ足してください。

STERI-CYCLE i160 LK 及び STERI-CYCLE i250 LK の水量:3|

運転を開始

- 1。ガラス扉またはガスシェードをあける。
- 2。 プレフィルター (2/図 5-1) を取り出す。
- 3。インサートプレート 注ぎ足しライン付き(STERI-CYCLE i160 LK): カバーを開けます (3/図 5-1)。
 非分割インサートプレート(STERI-CYCLE i250 LK): 容器の注ぎ足し作業に場所が足りない場合は、下のインサートプレート 取り出します。
- 4。充分に浄化した水を、装置内使用空間のタンク (蓋の下、ポジション 1 図 5-1) に 注ぐ。



図 5-1。 貯水槽

インキュベーターへの給水



図 5-2。 インキュベーターの排水・給水バルブ

インキュベーターの給水バルブから給水用漏斗で水を注ぐことができます。



図 5-3。 給水用漏斗

上図のようにガラスドアのロックに給水用装置のフックを引っ掛けます。 ホースを排水・給水バルブに接続します。

連通管の原理に従って、給水用漏斗によってインキュベーターの水位の高低差が決定されます。

給水用漏斗の水位線は、最小が0.5Lで、最大が3.5Lです。漏斗の最大容量は約0.4Lです。

- · インキュベーターの排水・給水バルブに漏斗を挿入します。
- 漏斗に少量の滅菌蒸留水を注いで、水位が安定するまで待ちます。これを数回繰り 返します。
- ・ 最大水位線に達するまで給水を繰り返します。

5。水量は、最大水量マーク「MAX」(1/図 5-4)を超えてはいけません。貯水槽(5/図 5-4)の最大水量は30です。



図 5-4。 MAX と書いてある注水マーク

- 6。多すぎる場合は、貯水槽の蓋で拭ってください。
- 7。プレフィルター(2/図 5-1)を再装着します。
- 8。下のインサートプレートを再装着します。もしくはカバーを閉じます。
- 9。CO₂/O₂/N₂供給設備のバルブが必ず開いているよう確実にしてください。
- 10。装置を電力スイッチでオンにする。
- 11。温度と CO₂/O₂ 含有量のの設定値を iCan ™タッチスクリーンで設定します。

装置スタート:

- 12。装置を自動スタートでスタートさせます(参照「自動スタートをオンにする」上 頁 6-14)。
- 13。経過表示 自動スタートがディスプレイに表示され、自動のスタートモードが始まり ます。
- 14。温度調整機が、温度設定値を調整し、湿度が発生します。
- 15。温度と相対湿度が一定になったら、CO₂/O₂計測システムとの整合が行われます。
- 16。CO₂/O₂ 調整は、設定した CO₂/O₂ 設定値に沿って行われます。
- 17。自動スタートモードが終わると、経過表示がオフになり、メインメニューに変わりま す。装置は運転可能です。

装置での処理:

18。装置内使用空間に培養菌をローディングします。

留意

自動スタートモードにかかる時間:

装置が冷たく周辺温度が低い場合、自動スタートモードに10時間かかる可能性があります。

ローディング:

サンプルにおける空気の循環と熱の均等を可能にするため、装置内使用空間のロー ディング面積は最大 70 % までの使用にとどめること。横に面積の大きい物や装置内 使用空間で熱排出装置を使用すると、熱の近郊に影響を及ぼす場合があります。横に 面積の大きい物や装置内使用空間で熱排出装置を使用すると、熱の近郊に影響を及ぼ す場合があります。

5 運転 インキュベーターへの給水



コンテンツ

- ・「電力スイッチ」上 頁 6-2
- ・「コントロールパネルと操作方法」上頁 6-3
- ・「iCan [™]タッチスクリーンの初期設定」上 頁 6-7
- ・「ループセンサーのウォームアップ」上頁 6-7
- ・「設定時のキーボタン操作」上 頁 6-8
- ・「温度設定値を設定する」上 頁 6-8
- ・「CO2 設定値を設定する」上 頁 6-9
- ・「02 設定値を設定する」上 頁 6-10
- ・「自動スタート機能」上 頁 6-12
- ・「Steri-Run を呼び出す」上 頁 6-16
- ・「ユーザーコンフィギュレーション」上頁 6-17
- ・「経過表示スケール」上 頁 6-43
- ・「エラー表示」上 頁 6-45
- ・「停電後の処置」上 頁 6-47

電力スイッチ



図 6-1。 電力スイッチ

電力スイッチは装置側面にあります。

- ・装置をオンにする:電力スイッチをおし、ライトが付きます。
 - ・短いシグナル音の後、ディスプレイが一時暗くなります。
 - ・ループのセンサーがウォームアップを始めます(「ループセンサーのウォーム アップ」上頁 6-7)。
- · 装置をオフにする: 電力スイッチおし、ライトが消えます。

コントロールパネルと操作方法

コントロールパネルはタッチスクリーン(iCan[™] タッチスクリーン)で、指やペンで軽 く触ることで操作ができます。



図 6-2。 メインディスプレイ: タッチディスプレイ

コントロールパネルの以下のディスプレイはタッチで操作が可能です

- ・ 温度表示 T,
- · CO₂表示 CO₂、
- ・ 表示 水位 RH,
- · 0₂表示(オプション)、
- 運転状況を示すアイコンシンボル、オプションのショートカットアイコン(参照 「アイコン説明」上 頁 6-40)、
- ・steri-run ボタン、
- ・メニューボタン、
- · 自動スタートボタン。

留意 タッチ操作域の拡大: エラー表示の確定をすると、タッチディスプレイ全体でタッチ操作が可能となります。

0₂/N₂ 調整なしの装備:



図 6-3。 iCan ™タッチスクリーン ガス供給 0₂/N_{2 なし}

コントロールパネルの機能ボタンと実測値表示、装置装備ガス供給 0₂/N₂ なしの場合:

- 1。表示部分 作業場温度実測値 (大きい数値が中央)、設定値 (小さい数字右上)と単位 (右下)
- 2。表示部分 CO₂ 濃度実測値 (大きい数値が中央)、設定値 (小さい数字右上) と単位 (右下)
- 3。HEPA フィルターの表示(その他のオプションシンボル参照 「アイコン説明」上 頁 6-40)
- 4。Steri-Run 除染モードスタートボタン
- 5。メニューナビゲーションを開けるボタン
- 6。自動スタートモードスタートボタン
- 7。表示 水位 RH

CO₂/O₂/N₂ マルチガス (オプション):



図 6-4。 iCan[™] タッチスクリーン、コンビガス接続付き

コントロールパネルの機能ボタンと実測値表示コンビガス供給 CO₂/O₂/N₂ 装備の場合:

- 1。表示部分 作業場温度 実測値 (大きい数値が中央)、設定値 (小さい数字右上) と 単位 (右下)
- 2。表示部分 CO₂ 濃度実測値 (大きい数値が中央)、設定値 (小さい数字右上)と単位 (右下)
- HEPA フィルターの表示(その他のオプションシンボル参照「アイコン説明」上 頁 6-40)
- 4。Steri-Run 除染モードスタートボタン
- 5。メニューナビゲーションを開けるボタン
- 6。自動スタートモードスタートボタン
- 7。表示部分 0₂ 濃度 実測値 (大きい数値が中央)、設定値 (小さい数字右上) と単位 (右下)
- 8。表示 水位 RH

操作方法

操作には、3段階あります:

- · A: ループ設定の直接操作: 温度、CO₂ 設定値、O₂ 設定値、
- · B: Steri-Run および自動スタートモードの開始、
- · C: 装置コンフィギュレーションにおけるサブメニューからのナビゲーション



図 6-5。 メニュー構造一覧

iCan [™]タッチスクリーンの初期設定

装置は、配送時には下記の設定値に設定されています:

- · 温度: 37°C
- · CO₂ 含有量: 5.0 %
- ・ 0₂ 含有量 (オプション): 21.0 %

留意 CO₂/O₂ 調整: 空気中の CO₂ 濃度が0%に近い値の場合、CO₂ 調整、およびループのエラーアラーム は設定値が0%の場合オフとなります。 空気中の O₂ 濃度が 21%のため、O₂ 調整およびループのエラーアラーム は、設定値 21%の際はオフとなります。

ループセンサーのウォームアップ

装置をオンにすると、ループセンサーが開始プロセス中にウォームアップに入ります。 ウォームアップの長さは様々です:

- 1。ループ 温度: 10 秒
- 2。ウォームアップ時間 CO₂ センサー (TCD と IR): 5分
- 3。0₂ループ:5分



図 6-6。 ウォームアップの表示

開始プロセスはシグナル音で通知されます。ウォームアップ中は、表示部分が数値の代わりに点(...)となります:

- · 温度表示、
- CO₂ 表示および
- · 0₂表示

ウォームアップ完了後、ループの実測値が表示されます。

留意 CO_2 ガス: O_2 ループの 5 分間にわたるウォームアップ後、装置内空間の CO_2 ガス放出および CO_2 ループの監視は行われません。

設定時のキーボタン操作

数値は、キーを押すことで段階的に高くしたり低くしたりすることができます:

・ - キーまたは + キー長押しをすることで、数値は高速で変化していき、

・約3秒間長押しを続けると、さらにその速度がアップします。

留意 設定の保存:

数値を変更後 Enter キーを押すと、その設定が保存されます。

設定を元に戻す:

30 秒間操作をしない状態 (タッチ操作が可能な部分およびキー)が続くと、メ ニューモードが自動的に 終了し、最後に保存した設定が有効となります。

温度設定値を設定する

1。温度キーを押します。
 ・温度メニュー(図 6-7)が表示されます。



図 6-7。 温度表示と温度選択メニュー

- 2。温度メニューを終了する:
- · 終了キー を押します。
- 3。温度設定値を設定する:
- · 設定値キーを押します。



図 6-8。 温度設定値を設定する

設定値 を上げる: ・+キーを押します。

設定値 を下げる:

- ーキーを押します。
- 4。設定値を保存する:
- · Enter キー を押します。
- ・メインメニューに戻ります。装置内空間にて計測された現在の実測値が、温度表示 部分に表示されます。

CO2 設定値を設定する

CO₂ 表示キーを押します。
 CO₂ メニューが表示されます。



図 6-9。 CO2 表示と CO2 メニュー

2。CO₂メニューを終了する: ・**終了**キーを押します。

Thermo Scientific

- 3。CO₂ 設定値を設定する:
 ・設定値キーを押します。
- 設定値 を上げる:
 ・+キーを押します。

設定値 を下げる: ・ ーキーを押します。

- 4。設定値を保存する:
- · Enter キー を押します。
- ・メインメニューに戻ります。装置内空間にて計測された実測値が、CO₂表示部分に表示されます。



図 6-10。CO₂ 設定値を設定する

留意 CO₂ループ オフ:

CO₂ 調整をオフにするためには、設定値を0%にします。 ループがオフになると、エラーアラームも作動しません。

ガスモニタリング:

CO₂ループの作動状況は、オプション機能のガスモニタリングには影響しません。CO₂ 調整がオンになっている場合でも、ガスモニターシステムはオンになっています。

留意

CO₂-設定値変更の後、内側容器を換気してください。これによって、アラームシステムの警告を避けることができます。

02 設定値を設定する

この設定は、オプションの $0_2/N_2$ 調整付き装備の場合のみ可能です。

1。ボタン 0₂ 表示を押します。

· 0₂ メニュー が表示されます。



図 6-11。02 表示箇所と02 メニュー

- 2。02メニューを終了します:
 ・終了キーを押します。
- 02 設定値を設定する:
 設定値キーを押します。

設定値 を上げる: ・+キーを押します。

設定値 を下げる:

- ・-キーを押します。
- 4。設定値を保存する:
- · Enter **キー** を押します。
- ・メインメニューに戻ります。装置内空間にて計測された実測値が、0₂表示部分に表示されます。



図 6-12。0₂ 設定値を設定する

留意

02-設定値変更の後、内側容器を換気してください。これによって、アラームシステムの警告を避けることができます。

留意初期設定: 0₂ センサーの型によって 0₂ 制御範囲のうち片側は工場で既に設定が行われています: 制御範囲 I:1 % - 21 % 制御範囲 II:5 % - 90 %

プロセスガスの使用:

 0_2 設定値が 21 % より低い場合、本装置を窒素供給に接続してください。 0_2 設定値が 21 % を超えている場合、本装置を酸素供給に接続してください。 $_0^2$ 設定値 21 % の場合、ループはオフになります。すなわち、エラーアラームについて も作動しません。ガスモニターシステム(オプション)は、それでもオンのままです。

留意

センサーがソケットに正しく差し込まれているかを確実にしてください。センサーが きちんと設置されていない場合、接触部に腐食が発生し自動スタートモード中に校正 値エラーが出る可能性があります。これは、センサーをオンにするだけで確認ができ ます。10 分経ってもエラー表示が現れない場合、自動スタートモードを開始してい ただけます。

自動スタート機能

自動スタート機能は、開始における自動化されたモードで、CO₂ 計測システムの調整を 行います。開始後、装置制御が温度を設定値に調整し、同時に湿度環境を作り出しま す。温度と相対湿度の値が安定したら、CO₂ 計測システムが自動でこの値への調整を行 い、設定の CO₂ 値で装置内使用空間ヘガスを送ります。 自動スタートモード使用上の注意:

CO₂ 計測システムの正確さを保つため、以下の場合は、本装置は常に自動スタートモードから開始してください:

- ・ 温度の設定値を1°C以上変更する場合、
- ・ 低湿度機能をオン / オフに変える場合、
- ・本装置が、ある程度の期間を置いた後に改めて使用される場合。

洗浄・メンテナンスの一環として、自動スタートモードは少なくとも四半期毎に運転 をしてください。

自動スタートモードにかかる時間:

自動スタートモードには、通常5~7時間かかります。室温が低い場合、そして装置 が温まっていない場合は、自動スタートモードに10時間かかることもあります。自動 スタートモード運転中にガラス扉を開けたり本装置への電力供給が遮断された場合、 このモードは中断されます。ガラス扉が再び閉められた場合、もしくは電力への再接 続が行われると、自動的に新たに開始されます。

自動スタートモード開始条件:

自動スタートモードの開始時は、装置内使用空間内の環境は周辺空気の値と同様です。 開始前に CO₂ および O₂ の設定値をご希望の値に設定しておいてください。装置内利用 空間の貯水槽には、充分な量の水が入っている必要があります。

自動スタートモード開始不能:

自動スタートモードは、以下のエラー発生時には開始することができません。

ループ 温度 :

- センサー破損、
- ・実測値が設定値を超えている(差異が大きすぎる)、
- ・実測値が設定値を下回っている(差異が大きすぎる)、
- ・実測値 妥当ではない
- ・ 校正値 大きすぎまたは小さすぎる
- ・ センサーへの接続が認識されない

ループ CO2 ガス供給:

センサーへの接続が認識されない。
 自動スタートボタンはこのようなエラー発生時には消え、押すことができません。

自動スタートモードのエラーによる中断:

自動スタートモードは、以下の場合中断されます:

- ・ 温度ループにおけるエラーが発覚した場合、
- CO₂ ループにおけるエラーが発覚した場合、
- 水の残量が少なすぎる場合、
- ・ 設定の CO₂ 値が許容範囲外の場合。

自動スタートモードを水なしで実行する:

自動スタートモードを水なしで、つまり装置内使用空間の貯水槽に水を入れずに実行 する場合、開始前に水センサーをオフにしてください(章「オプション」上 頁 6-31)。

自動スタートをオンにする

開始前の準備:

- 1。CO₂/O₂/N₂ガス供給設備のバルブが開いていることを確認します。
- 2。充分に浄化した水を、装置使用空間内の貯水槽に入れてください。注水量は、上の注 水マークを超えてはいけません。

自動スタートモードを呼び出す:

- 1。自動スタートボタンを押します。
- ・ 自動スタートプロセスのメニューが表示されます。



- 図 6-13。自動スタートをオンにする
- 2。自動スタートプロセスのメニューを終了し、自動スタートを中断:
- · 終了キー を押します。
- 3。必要に応じて水を足します。
- 4。自動スタートを開始:
- · Start ボタン を押します。

- 5。時間ブザーが鳴るまで 30 秒間、装置内使用空間を換気し、装置のドアを両方とも開 けます。
- 6。その後、装置のドアをどちらも閉めます。
- · 経過状況を示すステータスの表示が現れます。



図 6-14。自動スタートのステータス表示

```
留意中断:
自動スタートモードは、いつでも中断させることができます!
停止 ボタン を押します。
自動の再スタート:
自動スタートモードが以下の行動により中断された場合は、その後自動で再スタートされます:
・ ガラス扉を開ける、
・ オプションのガスシェードの場合は、外扉を開けた時、
・ 電力供給の中断
```

自動スタートを停止する

ステータス表示画面で**停止**ボタンを押すと、自動スタートモードを停止することができ ます。安全データ確認としてダイアログボックス自動スタート**停止**が表示されます。こ こで完全に停止させてしまうことも、再開することもできます。

- 1。自動スタートを再開する:
- · **戻る**キーを押します。
- · ステータス表示画面に戻り、自動スタートモードが再開されます。
- 2。自動スタートを停止する:
 - ·終了キーを押します。
 - ・

 中断の注意として三角の警告マークが、シグナル音と共に登場します。



図 6-15。自動スタートを停止する

- 3。中断の注意を承認する:
- ディスプレイを押します。どこでも構いません。
- · エラーダイアログボックス が出てエラー名が表示されます。

エラー システム - エラー: 自動スタート

図 6-16。自動スタート停止後のエラー表示

- 4。エラー表示を承認する:
- · 終了キー を押します。
- ・メインメニューに戻ります。

Ster i-Run を呼び出す

Steri-Run は、自動の除染モードで、装置内空間の消毒を行います。Steri-Run は、自動の除染モードで、装置内空間の消毒を行います。機能プロセスの詳細は、洗浄と消毒 についての章にてご確認ください(章8、ページ1)。

ユーザーコンフィギュレーション

ユーザーコンフィギュレーションを設定することにより、ユーザーインターフェースと 装置の追加機能を、日々の装置運転時のニーズに適合させることができます。ユーザー コンフィギュレーション選択メニュー(図 6-17)は、メインディスプレイにあるメ ニューボタンから呼び出すことができます。



図 6-17。ユーザーコンフィギュレーションメニュー

ユーザーコンフィギュレーションメニューは、6 つのカテゴリに分かれています:

- ・ 設定 / セットアップ,
- · イベント・ログ,
- ・オプション,
- ・アイコン説明,
- ・キーロック、
- ソフトウェアのバージョン。

ダイアログボックスにてユーザー独自の設定を行うには、図に示されているサブメ ニューに従い、ダイアログボックスを呼び出します。

設定 / セットアップ

選択メニュー 設定 / セットアップ(図 6-18)にて、入力ダイアログの数々にアクセス することができ、個々にあったユーザーインターフェースやポートのコンフィギュレー ションが設定できます:

- ・キーロックのコードを変更する。
- ・ 日付 / 時刻 を設定する,
- ディスプレイの明るさを設定する、
- ・キー音を設定する。
- ポートを設定する、
- ・ユーザーインターフェースの言語を設定する、
- リマインダーを設定する。



図 6-18。メニュー 設定 / セットアップ

キーロックのコードを変更する

キーロックを設定することで、権限のない者による設定の変更を防ぎます。この設定 で、数値の変更に関わるキーボタンのみがロックされます。 キーロックのコードは、4桁の数字です。

・発送時の初期設定は: 0000。

この設定は、ユーザー任意のコードに変更することができ、キーロックのダイアログ ボックス(「キーロックをオン/オフにする」上頁 6-42)にて変更をオンにすること ができます。

キーロックのコードを変更する:

1。**メニュー**キーを押します。

2。メニューコマンドのキーロックコードを選択する。 ·図 6-19 にある入力ダイアログが表示されます。

キーロックコード		
	コード:	
	1 2 3 4 5 6 7 8 9 0	
	削除 Enter 戻る	

図 6-19。キーロックのコードを変更する

- 3。初期設定の0000を入力:
 - ・ 新しく設定するコードの数字キーボタンを押し入力する。
- · 入力した数字コードは、隠れた状態で表示されます。

- 4。入力を確定する:
- · Enter キー を押します。
- · ディスプレイに、入力画面コード ニューと表示されます。空欄が4つあり、そこに 新しいキーコードが入力できるようになっています。

1 2 3 4 5 6 7 8 9 0		
Enter <<		

図 6-20。キーロックのコードを変更する

新しい4桁のコードを入力する:

- ·新しく設定するコードの数字キーボタンを押し入力する。
- · 数字コードが入力欄に表示されます。
- 5。数字を変更するために、カーソルを左に向けます: ・**戻す**(<<) キーを押します。
- 6。入力内容の適用し保存する:
- · Enter **キー** を押します。
- · 設定 / セットアップ選択メニューに戻ります。

(T

・メインメニューのアイコンバー(図 6-2 上 頁 3)に、キーロックが有効になっている印として、鍵マークが表示されます。

留意 一度設定したコードを変更する: 一度設定したコードは、同様の方法で変更することができます:

- · コード変更には、まずは現行のコードを入力し、
- 新しいコードを入力し確定します。

日付 / 時刻 を設定する

日付と時刻を、入力ダイアログで、ご希望のタイムゾーンに合わせて設定することができます。

- 1。**メニュー**キーを押します。
- 2。メニューコマンド 日付 / 時刻を選択します。
- 図 6-21 にある入力ダイアログが表示されます。

日付/時刻	
	日付
	時刻

図 6-21。選択メニュー 日付 / 時刻

3。日付の変更には、オプションの中から日付を選択します。 ・図 6-22 にある入力ダイアログが表示されます。



図 6-22。日付を設定する

- 4。日付を入力する:
- 数字キーを押します。
- 入力した数字が入力欄に表示されます。
- 5。数字を変更するために、カーソルを左に向けます: ・**戻す (<<)** キーを押します。
- 6。入力内容の適用し保存する:
- · Enter キー を押します。
- 7。日付 / 時刻の選択メニューに戻ります。
- 8。時刻の変更をするには、オプションの中から時刻を選択します。
- 図 6-23 にある入力ダイアログが表示されます。


図 6-23。時刻を設定する

- 9。時刻を入力する:
- ・ 数字キーを押します。
- 入力した数字が入力欄に表示されます。
- 10。数字を変更するために、カーソルを左に向けます: ・ **戻す** (<<) キーを押します。
- 11。入力内容の適用し保存する: • Enter キー を押します。
- 12。日付 / 時刻の選択メニューに戻ります。

ディスプレイの明るさを設定する

ディスプレイの明るさは、入力ダイアログにて1~100%の値の間で設定ができます。

明るさを設定する

- 1。**メニュー**キーを押します。
- 2。メニューコマンド 設定 / セットアップを選択します。
- 3。選択肢にあるディスプレイの明るさを選択します。
- ・ 🗵 6-24 にある入力ダイアログが表示されます。



図 6-24。ディスプレイの明るさを設定する

- 4。数値を上げる:
- ・ + キーを押します。
- 5。数値を下げる:
- ・-キーを押します。
- 6。値の変更は、表示欄に表示されます。ニューという表示になると、値に変更があった という意味ですが、設定の変更はまだ行われていない状態です。
- 7。変更の適用し保存する:
- Enter キー を押します。
- · 設定 / セットアップ選択メニューに戻ります。

キー音を設定する

キーボタンを押すときに鳴るキー音の音量は、入力ダイアログにて設定できます。 設定値は、0~100です。変更は5%ずつ可能です。

キー音の音量を設定する

- 1。**メニュー**キーを押します。
- 2。メニューコマンド 設定 / セットアップを選択します。
- 3。オプションの中からキー音を選択します。
- 図 6-25 にある入力ダイアログが表示されます。



図 6-25。キー音の音量を設定する

- 1。数値を上げる:
- ・ + キーを押します。
- 2。数値を下げる:
- ・ -キーを押します。
- 3。値の変更は、表示欄に表示されます。ニューという表示になると、値に変更があった という意味ですが、設定の変更はまだ行われていない状態です。
- 4。変更の適用し保存する:
- Enter キー を押します。
- · 設定 / セットアップ選択メニューに戻ります。

USB インターフェースのボーレートを設定する

USB インターフェースのデータ移行率は、入力ダイアログにて設定できます:

インターフェースのデータ移行率は、既定のボーレート(9600、19200、38400、57600 ボー)内で設定を変更することができます。

ボーレートを設定する

- 1。**メニュー**キーを押します。
- 2。メニューコマンド 設定 / セットアップを選択します。
 - · 図 6-26 にある入力ダイアログが表示されます。



- 図 6-26。選択メニュー インターフェース USB
- 3。オプションの中からインターフェース USB を選択します(図 6-26)。 ・図 6-27 にある入力ダイアログが表示されます。



図 6-27。USB インターフェースのボーレートを設定する

- 4。+またはーキーで、USB インターフェースのボーレートを合わせます(図 6-27)。
- ・ボーレートを上げる:+キーを押します。
- ボーレートを下げる: ーキーを押します。
- 5。値の変更は、表示欄に表示されます。ニューという表示になると、値に変更があった という意味ですが、設定の変更はまだ行われていない状態です。
- 6。変更の適用し保存する:
- Enter キー を押します。
- · 選択メニュー インターフェース USB に戻ります。
- 7。新しい設定を有効にする:
- ・メインメニューに戻ります。
- · 約 10 秒待ち、再起動し、装置を電力スイッチでオフ / オンにします。

ユーザーインターフェースの言語を設定する

ユーザーインターフェースの言語を入力ダイアログで設定します。7 つの言語からお選びいただけます:

- ・ ドイツ語、
- ・英語、
- ・スペイン語、
- フランス語、
- ・イタリア語、
- ・中国語、
- ・日本語。

表示言語を選択する:

- 1。メニューキーを押します。
- 2。メニューコマンドで言語を選択します。
- 図 6-28 にある入力ダイアログが表示されます。



図 6-28。言語を設定する

- 3。選択肢の中で上にページをめくる:
- ・ + キーを押します。
- 4。選択肢の中で下にページをめくる:
- ・ -キーを押します。
- 5。新たに選んだ言語が、表示欄に表示されます。ニューという表示になると、値に変更 があったという意味ですが、設定の変更はまだ行われていない状態です。
- 6。選択内容を適用し保存する:
- · Enter **キー** を押します。
- ・ 設定 / セットアップ選択メニューに戻ります。

リマインダーを設定する

リマインダーは、装置制御の通知・コントロールシステムの一部です。装置の二大重要 機能である Steri-Run および自動スタート、ならびに定期的なサービスのために、一定 の時期が来たら通知が行われるように設定できます。 時期の開始は、その前のリマインダー発動した日の 00:00 時からです。

リマインダー発動すると、ディスプレイに通知メッセージが表示されます:

- Steri-run: Steri-run を実行してください。
- ・ 自動スタート: 自動スタートを実行してください。Steri-Run 除染モードが完了したら表示されます。
- ・サービスインターバル:サービスを実施してください。サービス通知を確定してく ださい。アイコンサービス請求が以下に表示されます。

通知メッセージは、それぞれの事項が実施完了したら、消えます。

初期設定

90 日
オフ
Aオフ
365日

リマインダーを設定する

- 1。**メニュー**キーを押します。
- 2。メニューコマンド リマインダーを選択します。 ・図 6-29 にある入力ダイアログが表示されます。



- 図 6-29。リマインダー機能を選択します。
- 3。該当のメニューコマンド、例えば steri-run を選択します。
- ・ 図 6-29 にある入力ダイアログが表示されます。



図 6-30。steri-run のインターバルを設定する

- 1。日にちを増やす:
- ・+キーを押します。
- 2。日にちを減らす:
- ーキーを押します。
- 3。値の変更は、表示欄に表示されます。ニューという表示になると、値に変更があった という意味ですが、設定の変更はまだ行われていない状態です。
- 4。インターバルをオフにする:
- ・数値を**オフ**に合わせる。
- ・ -キーを押します。
- 5。変更の適用し保存する:
- · Enter キー を押します。
- ・選択メニューリマインダーへ戻ります。

イベント・ログ

選択メニュー データ記録 (図 6-31) では、装置運転中のイベントの記録およびアウト プットのダイアログにアクセスすることができます:

- ・イベント表示,
- ・インターバル(記録インターバル)、
- ・エラー表。



図 6-31。選択メニュー データ記録

イベントを表示する

イベント表示は、装置運転中に記録されたイベントが、日付と時刻と共に、端的に1つ 1行の形での報告です。

情報は、時間順に、最新のものが上に来る形のリストとなっています。リストは見ることはできますが、編集はできません。イベント表示がすうページに及ぶ場合は、ディス プレイ内でページをめくることができます。ステータス表示で、現在何ページ目にいる かを知ることができます。

イベント表示を呼び出す

- 1。**メニュー**キーを押します。
- 2。メニューコマンド データ記録を選択します。 ・図 6-31 にある入力ダイアログが表示されます。
- 3。メニューコマンド イベント表示を選択します。
 - ・ 🗵 6-32 のリストが表示されます。



図 6-32。イベントを表示する

- 1。リスト内でページを進める:
- ・**次**キーを押します。
- 2。リスト内で前のページに戻る:
 - ・ **前**キーを押します。
- 3。表示を終了する:
- · **終了**キー を押します。
- ・選択メニュー データ記録 に戻ります。

記録インターバルを設定する:

保存容量が決まっているため、新しいイベントが記録されると、一番古い物が削除され ます。記録されたイベントの時期は、ご選択の記録インターバル時間に大きく関わって きます。

イベント表示を呼び出す

- 1。**メニュー**キーを押します。
- 2。メニューコマンド データ記録を選択します。 · 図 6-31 にある入力ダイアログが表示されます。
- 3。メニューコマンド 記録インターバル時間を選択します。 ・図 6-33 にある入力ダイアログが表示されます。



図 6-33。記録インターバルを設定する

記録インターバル	最大表示時間
10 秒	22.5時間
30 秒	2.8 日
60 秒	5.6 日
120 秒	11.2 日
180 秒	16.8 日
3600 秒	336 日

装置運転中のループ計測値を記録し、経過の再表示がされる記録インターバルは、秒単位で設定することが可能です(「経過表示スケール」上頁 6-43)。

10 秒~3600 秒の範囲内で設定が可能です。

- 1。数値を上げる:
- ・ + キーを押します。
- 2。数値を下げる:
- ーキーを押します。
- 3。値の変更は、表示欄に表示されます。ニューという表示になると、値に変更があった という意味ですが、設定の変更はまだ行われていない状態です。
- 4。変更の適用し保存する:
- Enter キー を押します。
- · 選択メニュー データ記録に戻ります。

留意データ記録の保存サイクル:

記録インターバル時間は、エラー表の記録に影響は及ぼしません。

エラー表を見る

エラー表は、装置内のコントロールシステムが発見したエラーを、最新のものから順に リスト形式で表示したものです。一番最近検出されたエラーが、一番上に表示され、全 部で22つの記録があります。記録一つには、エラーが発見されたループ、日付、時刻 およびエラー名が含まれています。エラー表は見ることはできますが、編集はできませ ん。イベント表示が2ページにわたる場合、ディスプレイ内でページをめくることがで きます。ステータス表示 001/002 または 002/002 は、現在のページを示しています。

エラー表を見る

- 1。**メニュー**キーを押します。
- 2。メニューコマンド データ記録を選択します。
- 図 6-31 にある入力ダイアログが表示されます。
- 3。メニューコマンド イベント表示を選択します。
- ・ 図 6-34 のリストが表示されます。



図 6-34。エラー表を見る

留意 エラー対応: エラーの原因やエラー解消についての詳細一覧は、本章の終わりにあります!

- 1。エラー表内でページを進める:
- ・ **次**キーを押します。
- 2。リスト内で前のページに戻る:
- ・ **前**キーを押します。
- 3。表示を終了する:
- · **終了**キー を押します。
- ・選択メニュー データ記録に戻ります。

オプション

選択メニュー オプション (図 6-35) 装置オプション機能の設定ダイアログ全てにアク セスすることができます:

- ・アラーム、
- ・低湿度、
- ガスシェード (オプション)、
- ・ 空気湿度センサー、
- ・ブザー音,
- ・0₂ガス供給(オプション)、
- ・HEPA フィルター、



図 6-35。選択メニュー オプション

アラームリレーを設定する

アラームリレーは、装置内部コントロールシステムを外部の電気供給アラームモニタリ ングシステムにつなぐ際のインターフェースとなります。外部モニタリングシステムの 入力シグナルに応じて、電力供給アラームはオンまたはオフにすることができます。電 力供給アラームがオンの場合、停電はエラーとして認識されます。アラームリレーのワ イヤリングについては「アラームコンタクトを接続:」上 頁 4-17 章に記載があります。

- 1。メニューキーを押します。
- 2。メニューコマンド オプションを選択します。
- · 図 6-35 にある入力ダイアログが表示されます。
- 3。メニューコマンド アラームを選択します。
- ・図 6-36 にある入力ダイアログが表示されます。



図 6-36。選択メニュー アラーム

- 4。メニューコマンド アラームリレーを選択します。
- · 図 6-37 にある入力ダイアログが表示されます。

アラームリレー			
	電力供給ア	' ラーム	
	現在	オン	
	設定	オン	
	-	Enter	+

図 6-37。アラームリレーを設定する

- 1。二つのコンディションの間で切替える:
- ・ + キーを押します。

または

- ーキーを押します。
- 2。変更の適用し保存する:
- · Enter キー を押します。
- ·選択メニューオプションに戻ります。

低湿度を設定する

相対湿度が高いために培養容器にコンデンゼーションが形成されてしまう場合、装置内 使用空間の湿度を下げることができます。発送時は、装置操作は高湿度、つまり相対湿 度約 93 %に設定されています。

低湿度をオンにしている状態では、装置内使用空間の相対湿度は約 93 % から約 90 % に 下がります。

変更をする場合は、調整に時間がかかる可能性があります。培養容器のコンデンゼー ションを効果的に避けるために、継続的にこの設定にするようにしてください。

装置内空間の空気湿度を下げる

- 1。**メニュー**キーを押します。
- 2。メニューコマンド **オプション**を選択します。
- ・図 6-35 にある入力ダイアログが表示されます。
- 3。メニューコマンド 低湿度を選ぶ。
- 図 6-38 にある入力ダイアログが表示されます。

低湿度			
	低湿度		
	現在	オフ	
	設定	オフ	
	_	Enter	+

図 6-38。低湿度を設定する

- 1。二つのコンディションの間で切替える:
- ・+キーを押します。 または
- ・-キーを押します。
- 2。変更の適用し保存する:
- · Enter キー を押します。

選択メニュー オプションに戻ります。

70

メインメニューに戻ると、低湿度アイコンが表示されているのが確認できます。

留意低湿度: 低湿度機能をオン/オフにすると、イベントリストに記録されます。

ガスシェードを設定する

サンプルにアクセスする際の開口面が小さく作ってあるので、オプションのガスシェードが 搭載されている装置は、インキュベーションパラメータのインターバルも短く済みます:

- · 装置内使用空間温度、
- ・ CO₂ 濃度,
- ・0₂ 濃度、
- ・相対湿度。

装置装備変更をしたら、装置制御をオプション内ガスシェードに切り替える必要があります。

留意 誤作動: ガスシェードへの切り替えは、コントロールパラメータの変更に影響します。 ガスシェード機能が実際に使用されている扉の種類以外の物で設定されている場合、 インキュベーションに誤作動が発生する可能性があります。

ガスシェードを設定する

- 1。**メニュー**キーを押します。
- 2。メニューコマンド **オプション**を選択します。
- · 図 6-35 にある入力ダイアログが表示されます。

3。メニューコマンド ガスシェード を選ぶ。
 · 図 6-39 にある入力ダイアログが表示されます。

ガスブラインド			
	ガスブラ	インド	
	現在	ガラスドア	
	設定	ガラスドア	
	_		+

図 6-39。ガスシェードを設定する

- 1。二つのオプションの間で切替える:
- ・+キーを押します。
- または
- ・ -キーを押します。
- 2。変更の適用し保存する:
- · Enter キー を押します。
- ·選択メニューオプションに戻ります。

水位センサーをオン / オフにする

周囲湿度でのインキュベーション運転時や、自動スタートモードを水なしで行う場合、 水位センサーはオフにすることができます。それにより、装置内部コントロールシステ ムの水位センサー通知が行われなくなります:



注意 水位センサーがオフの時は、Steri-run モードが水でも開始することができてしまいますが、これは誤った使用でありファンモーターが故障する可能性があります。

- 1。メニューキーを押します。
- 2。メニューコマンド **オプション**を選択します。
- · 図 6-35 にある入力ダイアログが表示されます。
- 3。メニューコマンド水位センサーを選ぶ。
 - 図 6-40 にある入力ダイアログが表示されます。

水位センサー			
	水位セン!	ナー	
	現在	オン	
	設定	オン	
		Enter	+

図 6-40。水位センサーを設定する

- 1。二つのコンディションの間で切替える:
- · + キーを押します。
 - または
- ・ ーキーを押します。
- 2。変更の適用し保存する:
- · Enter キー を押します。
- ·選択メニューオプションに戻ります。

ブザー音 オン / オフにする:

装置内コントロールシステムがエラーを認識すると:

- エラーが表示されアラームリレーをオンになる他、
- ・ 聴覚アラームとしてブザー音が鳴る。

ブザー音は、ミュート設定ができます。

- 1。**メニュー**キーを押します。
- 2。メニューコマンド**オプション**を選択します。
- · 図 6-35 にある入力ダイアログが表示されます。
- 3。メニューコマンド ブザー音を選ぶ。
- · 図 6-41 にある入力ダイアログが表示されます。



図 6-41。アラームリレーを設定する

- 1。二つのコンディションの間で切替える:
- ・+キーを押します。

または

- ーキーを押します。
- 2。変更の適用し保存する:
- · Enter キー を押します。
- · 選択メニュー オプションに戻ります。

02調整をオン / オフにする

作業プロセスへの要求事項に応じて、02調整をオンまたはオフにすることができます。 この設定は、オプションの02/N2調整付き装備の場合のみ可能です。

- 1。**メニュー**キーを押します。
- 2。メニューコマンドオプションを選択します。
 図 6-35 にある入力ダイアログが表示されます。
- 3。メニューコマンド 02 を選ぶ。
 · 図 6-42 にある入力ダイアログが表示されます。



図 6-42。02 調整をオン / オフにする

- 1。0₂調整の二つコンディションの間で切替える:
- ・+キーを押します。 または
- ーキーを押します。
- 2。値の変更は、表示欄に表示されます。ニューという表示になると、値に変更があった という意味ですが、設定の変更はまだ行われていない状態です。
- 3。設定を適用し保存する:
- · Enter キー を押します。
- ·選択メニューオプションに戻ります。

留意 0₂ 値の表示: 0₂ 調整がオフの場合、0₂ 表示には実測値が表示されません(---)。 これにより、0₂ センサーの酷使を防ぎます。設定値が21%に設定されている場合、0₂ ループのアラームは作動しません。これは、0₂ 制御範囲両方に適応します: ・制御範囲 I:1%-21%

· 制御範囲 II: 5 % - 90 %

それから 02 表示に実測値が表示されます。

装置内空間を換気する

本装置が 0_2 または N_2 で運転される場合、 0_2 調整がオフになった後に装置内空間を換気 する必要があります。

ガスモニタリング

 0_2 ループのオンオフは、オプションのガスモニタリング機能に影響はありません。 0_2 調整がオフの場合や、 0_2 センサーが無効の場合も、ガスモニターシステムはオンになっています。

HEPA フィルターをオン / オフにする:

本装置にHEPA フィルターの取付けなしで運転を行う場合、誤作動防止のためコンフィ ギュレーションでオフ設定をする必要があります。

- 1。メニューキーを押します。
- 2。メニューコマンド **オプション**を選択します。
- ・ 🗵 6-35 にある入力ダイアログが表示されます。
- 3。メニューコマンド HEPA フィルターを選ぶ。
- ・ 図 6-43 にある入力ダイアログが表示されます。



図 6-43。HEPA- コンフィギュレーション

4。メニューコマンド HEPA コンフィギュレーションを選ぶ。 ·図 6-44 にある入力ダイアログが表示されます。



図 6-44。HEPA フィルターをオン / オフにする

- 1。二つのコンディションの間で切替える:
- ・+ キーを押します。 または
- ・-キーを押します。

- 2。変更の適用し保存する:
- · Enter キー を押します。
- ·選択メニューオプションに戻ります。

ISO ・メインメニューのアイコンバーに(図 6-2 上 頁 3)5分後に HEPA フィルターのオ 5 ン表示 ISO 5 が出ます。

アイコン説明

重要な運転状況やエラー表示、例えばキーロック、低湿度などは、データ記録やエラー 表に記録されるだけでなく、タッチディスプレイのメインメニューにアイコンとして表 示されます。アイコンの意味は、ダイアログボックス アイコン説明に説明があります。

アイコン説明を呼び出す

- 1。**メニュー**キーを押します。
- 2。メニューコマンド アイコン説明 を選ぶ。 · 図 6-45 にある情報ダイアログが表示されます。

アイコン説明				
	Ű-	ガスボトル 空	ISO	HEPA稼動
	ŀ	過温度	X	サービス 請求
	€	キーロック		
	7	低湿度		

図 6-45。アイコン説明

- 表示を終了する:
- · 終了キー を押します。
- · 選択メニュー ユーザーコンフィギュレーションに戻ります。

それぞれのアイコンの意味:

ガスボトル 空:

Ű.

◄

1 本もしくは複数のガスボトルの残量が少なすぎる旨を示すエラー表示で、規定通りの ガス供給を行うよう促すものです。 このアラーム機能は、ガスモニタリングシステム(オプション) 搭載の装置にのみあり ます。

過温度:

装置調整が過温度保護機能をオンにし緊急調整に切り替わったことを示すエラー表示です。

キーロック:



キーロックがオンにされ、設定変更ができない状態にあることを示す機能表示です(説明書参照「キーロックをオン/オフにする」上頁 6-42)。

低湿度:



装置内空間の相対湿度が約93%から約90%に下がったことを示す機能表示です(説明 書参照 「低湿度を設定する」上頁 6-33)。

HEPA フィルターオン:

ISO 5 装置内空間で HEPA フィルターがオンになったことを示す機能表示です(説明書 参照 「HEPA フィルターをオン/オフにする:」上 頁 6-39)。

サービス請求:



定期的なサービスリマインダーの時期が来たことをお知らせする通知です。アイコンの 表示は、リマインド通知が確認されると、ダイアログボックス リマインダーにて時期を 入力することで操作ができます。

ガスボトル残量の表示 (オプション):

本装置にガスモニタリング機能(オプション)が搭載されている場合、ガスボトルAと ガスボトルBのアイコンが、それぞれ CO₂ または O₂ メニュー に表示されます。アイコ ンはガスボトルの残量(満/空)を示します。青枠で囲われたアイコンは、次の供給 にどのガスボトルに切り替えができるかを示しています。



図 6-46。ガスモニタリングのアイコン

- ・ 満タンのガスボトルは、手動で交換できます。
- ・ 青枠で囲われたボトルのアイコンを押します または
- この場合は自動で実行されます:
- ・ボトル圧が 0.6 bar 以下に落ちた場合。

手動または自動でガス供給が切り替えられると、30秒の間は更なる切り替えができません。その後だいたい2分ほどで、交換ボトルの残量を新たに算出し、表示します。 ガスモニタリングシステムは、接続されているボトル両方の残量を管理しています。 ボトルが空になると:

- アラーム音は鳴らず、
- エラーが表示され、
- イベントリストに記録がされます。

ボトルが両方とも空の場合:

- アラーム音がなり、アラームリレーがオンになり、
- エラーが表示され、
- イベントリストに記録がされます。
- イベントリストに記録がされます。

留意 ボトルの交換:

自動もしくは手動のボトル切替えは、イベントリストに記録がされます。

キーロックをオン / オフにする

入力ダイアログで、キーロックのオンオフ操作が可能です。発送時は、キーロックは標 準コード 0000 に設定されています。

- 1。数字キーを使用し、4桁の数字を入力します。入力内容は、表示欄に暗号化し表示されます。
- 2。間違った入力を全て削除する:
 ・削除キーを押します。
- 3。入力を中断する:
- · **戻る**キーを押します。
- · 選択メニュー ユーザーコンフィギュレーションに戻ります。
- 4。入力を確定する:
- · 終了キー を押します。
- · 選択メニュー ユーザーコンフィギュレーションに戻ります。

留意 コードを変更する: 現行のコードを、ダイアログボックス キーロック / メニューコード 設定 / セット アップにて設定し直すことができます(「キーロックのコードを変更する」上 頁 6-18)。

コードのリセット: キーロックコードの標準コードへのリセットは、Thermo Fisher Scientificのテク ニカルサポートのみ行うことができます。



図 6-47。キーロックをオン / オフにする

ソフトウェアのバージョン

表示欄に、装置制御に使用されているソフトウェアのバージョンがひょうじされます。

・表示を終了する: ・ 終了 キー を押します。 ・選択メニュー ユーザー=	ュンフィギュレーション に戻り	ります。
バージョン		
	ソフトウェア メインボード: ソフトウェア CO2センサー: ld RHセンサー: ソフトウェア ディスプレイ: コンフィギュレーション:	50134596-0030 50138673-0001 50141134-0001 50130789-0026 8828454

図 6-48。ソフトウェアのバージョン

経過表示スケール

ループ3つの経過表示:

・温度,

- 0... 20% CO₂,
- 0...100% 0₂,

表示スケールは2パターンあります。

1。全画面表示

- · メインメニューの CO2 表示 ボタンを押します。
- · CO₂(図 6-9)メニューが表示されます。



図 6-49。CO2 濃度の経過表示 を呼び出す

2。アイコン 推移を表示 を押します。 ・経過の表示がされます。



図 6-50。CO2 濃度の経過表示 を呼び出す

3。拡大表示:

- ダイアグラムのご希望箇所を、指 / ペンで真四角を引きます。四角の大きさは、スタートポイント(プッシュポイント、ダイアグラム上枠左)を引っ張り、エンドポイント(放す、ダイアグラム下枠右)で定めます。
- ・ 選択した四角の中を、どこでもいいので押します。選択箇所が拡大表示されます。
- この動作は、拡大された箇所がご希望の大きさになるまで、または最大拡大レベル (最大 30 データログ分、記録インターバル 60 秒設定の場合は 30 分に相当)に到達 するまで、何度でも繰り返し行っていただけます。
- · ズームモード内で、ページをめくることもできます。

- 4。再び全体表示にする:
- ・ダイアグラムの小さい範囲の上絵にある正方形を引っ張り、マーキングされている ところ以外の画面をどこでもよいのでを押します。
- 5。経過表示を閉じる:
- · 終了キー を押します。
- ・メインメニューに戻ります。

留意 記録インターバル:

データ記録のインターバルは、ダイアログボックス 記録インターバル時間(「記録 インターバルを設定する:」上 頁 6-29)にて設定し直すことが可能です。

エラー表示

エラー検知は、装置内コントロールシステムの一部です。ループおよびそのセンサーの 監視を行います。システム内でエラーが報告されると、アラームリレーがオンになり、 以下のシグナルおよび通知がされます:

- ・聴覚アラームとしてブザー音が鳴る、
- 三角の警告マークが点滅し、該当のアイコンがメインメニューで表示される;そう すると値の表示が更新されない、
- 検知したエラーはエラー表に記録される、
- 経過はイベント表示に記録される。
- エラー確定後も尚エラーが消えない場合は、赤いボタン(T、CO₂、O₂、RH またはシ ステム、メニューキー)にて除くことができます。

エラー表示イベントに対するリアクション

アラームリレーが操作アクションによってオンになってしまった場合、エラー表示を確 定することで元に戻すことができます(例えば手動で Steri-Run 除染モードを中断した 時)。

アラームリレーがテクニカルエラーによってオンになると、エラーの原因が解消される までオンのままとなります(例えば、装置内空間の水位が低すぎるなど)。

- 1。エラー表示を承認する:
- · 三角の警告マークが表示されたら、タッチディスプレイに触れます。
- ・ダイアログボックス エラー が現れ、検知されたエラー原因が表示されます。
- ・シグナル音はオフになります。
- 2。エラー表示を閉じます:
 - · 終了キー を押します。
- エラー表示は消えます。



図 6-51。イベントエラー表示

過温度保護機能リセット



装置制御が過温度保護機能をオンにし緊急調整に切り替わると、三角の警告マークが点 滅しメインメニューにアイコンが表示されます。



図 6-52。エラー表示 過温度

- 1。エラーの原因を表示する:
- タッチディスプレイに触れます。
- ・ダイアログボックス エラー が現れ、検知されたエラー原因が表示されます。
- ・シグナル音はオフになります。

エラー 温度実測値が高すぎます	

図 6-53。エラー表示 過温度

- 2。エラー表示を閉じます:
- · 終了キー を押します。
- エラー表示は消えます。
- ・ 温度表示が赤で囲われます。
- 3。エラー表示を戻す: ・装置をオフにする。
- 4。扉を開けて、装置内空間の冷却をします。
- 5。装置をオンにする。

エラーの原因(参照 エラー表)を取り除いても尚、過温度保護機能がオンになってし まう場合は、本装置の運転を停止し、テクニカルサービスの請求を行ってください。

停電後の処置

電力供給中断の後、内部湿度のコンデンセーションがセンサー部分で発生する可能性が あります。これにより、センサーの機能が妨げられ、誤った測定値が表示されたり装置 が故障サインを出したりということが起こる可能性があります(センサー破損、「エ ラーの原因とエラー対処の一覧」上 頁 6-48 参照)。

誤りなく装置の運転ができるようにするために、以下の処置を行ってください。

- 1。排水し内部を乾かします。
- 2。装置を1時間水なしで 55°C に温めます。
- 3。その後装置の扉を開けて冷まします。
- 4。「装置使用の開始」上 頁 4-1 章の通りに、インキュベーションの温度で運転します。

他の方法としては、もしくは上記の処置が効かない場合は、180 °Cの消毒をスタートさせてください。これに関しては、章「Steri-Run 除染モード」上 頁 8-6 を参照ください。

約1時間後、消毒モードを中断することが可能です。これで、センサーは乾いたはずで す。これで、センサーは乾いたはずです。

エラーの原因とエラー対処の一覧

エラー表には、エラー源、エラー原因 およびエラー解消方法についての情報が記載し てあります。

Thermo Fisher Scientific テクニカルサービスとのやり取りのため、装置データを準備 しておいてください。

ルー プ	エラー表示	原因	処置	アラーム - リレー	ブザー音	ログ
システム	装置ドアの開扉時 間が長すぎ	装置ドアが 10 分以 上あいたまま	装置ドアを閉じる	Х	Х	Х
	エラー: ディスプ レイ	ディスプレイがメイ ンボードと反応しな い ^{*1)}	装置リセット。何 度も起こるようで あれば、サービス を請求する	X	Х	Х
	エラー:EEPROM メ インボード	メインボードの EEPROM に欠陥	装置リセット。何 度も起こるようで あれば、サービス を請求する	Х	Х	Х
	エラー:データログ	データログ保存の際 のエラー。インキュ ベータは、引き続き 機能可能です。	装置リセット。何 度も起こるようで あれば、サービス を請求する			
	エラー: steri-run	Steri-Run モード中 のエラー	装置リセット。何 度も起こるようで あれば、サービス を請求する	Х	Х	Х
	steri-run 中のパ ワーダウン	Steri-Run モード時 の停電	装置を再起動し、 Steri-Run を新た に開始する。	Х	Х	Х
	エラー: 自動ス タート	自動スタートモード 中のエラー	自動スタートを新 たに開始する。何 度も起こるようで あれば、サービス を請求する。	X	X	Х
	エラー: ADC	参考抵抗値が公差範 囲外	装置リセット。何 度も起こるようで あれば、サービス を請求する。	X	Х	Х
	エラー:ファン	ファンの実測値が公 差範囲外です。	装置リセット。何 度も起こるようで あれば、サービス を請求する。	Х	Х	Х
	IR センサー交換	新しいシリーズ番号 認識	アラーム停止	Х	Х	X

ルー プ	エラー表示	原因	処置	アラーム - リレー	ブザー音	ログ
温度	センサー破損	計測値が、許容リ ミット外	サービス請求。 センサーを熱して 乾かす。	Х	Х	Х
	実測値 大きい	実測値 > 設定値 + 1°C *2) *4)	許容 周囲温度を 超えない / サー ビスを請求。	Х	Х	Х
	実測値 小さい	実測値 < 設定値 - 1°C ^{*3)*4)}	エラーが自動的に 消えない場合は、 サービスを請求。	Х	Х	Х
	実測値 妥当ではな い	温度シグナルの妥当 性がない	装置リセット。何 度も起こるようで あれば、サービス を請求する。	Х	Х	Х
	校正値 大き / 小さ すぎ	温度の最大整合値を 超えた / 下回った	アラームを停止 し、ターゲット値 を入力する。			Х

ルー プ	エラー表示	原因	処置	アラーム - リレー	ブザー音	ログ
020% CO ₂	センサー破損	計測値が、許容リ ミット外	自動スタートを実 行する。その後引 き続きエラーが出 た場合は、「停電 後の処置」上 頁 6-47 章のエ ラー解決を参照。 エラーが解決しな い場合は、サービ スに連絡。	X	X	X
	実測値 大きい	実測値 > 設定値 + 1% ^{*4)}	自動	Х	Х	Х
	実測値 小さい	実測値 < 設定値 - 1% ^{*3) *4)}	自動	Х	Х	Х
	RH 接続障害	RH センサーがメイ ンボードと反応しな い	自動	Х	Х	Х
	校正値 大き / 小さ すぎ	CO ₂ の最大整合値を 超えた / 下回った	アラーム停止			Х
	接続障害	センサーがメイン ボードと反応しない	自動	Х	Х	Х
	エラー: ガスボト ル切替え装置	ガスボトル切替え装 置がメインボードと 反応しない	自動	Х	Х	Х
	ガスの不足	CO ₂ ガスボトルは両 方とも空です	少なくとも CO ₂ ボ トル 1 本交換す る。	Х	Х	Х
	RH センサー破損	計測値が、許容リ ミット外	サービス請求。 「停電後の処置」 上 頁 6-47 章も 参照。	Х	Х	Х
	ガスボトル A 空	ガスボトル A は空	ガスボトル A 交換			
	ガスボトル B 空	ガスボトル B は空	ガスボトル B 交換			

ルー プ	エラー表示	原因	処置	アラーム - リ レー	ブザー音	ログ
0100% 02	センサー破損	計測値が、許容リ ミット外	サービス請求	Х	Х	Х
	実測値 大きい	実測値 > 設定値 + 1% ^{*4)}	ガス供給を確認す る。最大 1 bar まで下げる。	Х	Х	Х
	実測値 小さい	実測値 < 設定値 - 1% ^{*4)}	ガス供給を確認す る。ガスボトルを 交換する 最大 1 bar まで上げる。 供給管を確認する。	X	X	X
	接続障害	センサーがメイン ボードと反応しない	サービス請求	Х	Х	Х
	エラー: ガスボト ル切替え装置	ガスボトル切替え装 置がメインボードと 反応しない	自動	Х	Х	Х
	ガスの不足	0 ₂ ガスボトルは両 方とも空です	少なくとも 0 ₂ ボ トル 1 本交換す る。	Х	Х	Х
	ガスボトル A 空	ガスボトル A は空	ガスボトル A 交 換			
	ガスボトル B 空	ガスボトル B は空	ガスボトル B 交 換			
rН	水なし	貯水槽内の水量が足りない。	水を注ぎ足すか、 求なしてする でする。水なしてす る。大変にす る。大変にす の た で た の い 場 合 に す て た な た た た た た た た た て に す の し て す の し て す の し て す の の の の の の の の の の の の の の の の の の	X	X	X

*1) エラーはディスプレイにのみ表示され、エラー表には記録されません。

*²⁾ エラーが発生したら、サンプル保護のために特別調整 がオンになります。これを明確にするため、アイコンも表示されます。

*3) エラー表示までの待ち時間:

・扉を開けてから45分

・設定値変更から 159 分

*4) この値は、サービス員により変更可能です。

<mark>6 運転</mark> 停電後の処置

装置の利用停止

コンテンツ

・「装置の利用を停止する」上 頁 7-1

装置の利用を停止する



注意 汚染の危険性があります! 装置内空間の表面は、汚染されている可能性があります。周囲に菌が移ってしまう危 険があります。 利用停止の際は、装置を除染すること!

- 1。培養容器と補助器具などを装置内空間から取り出す。
- 2。充分に容量のある回収容器を用意する。
- 3。ホースの先を回収容器に入れてホースバルブを安全に留める。
- 4。貯水槽が空になっていく。



- 図 7-1。 給水・排水 貯水槽の排水バルブ
- 5。貯水槽の水を全て回収容器に排出し、空にする。
- 6。ホースバルブを再び外す。
- 7。Steri-Run 除染モードを開始する(「Steri-Run を呼び出す」上 頁 6-16)。

- 8。Steri-run 除染モードが完了したら、装置を電力スイッチで切る。
- 9。電力プラグをコンセントの差込口から抜き、再接続を防ぐ。
- 10。CO₂/O₂/N₂ 供給設備のクローズバルブを閉じる。
- 11。ガス耐圧ホースを装置背面の接続バルブから抜く。
- 12。装置デコミッショニング中は装置内空間を常に換気状態にしてください。それにはガ ラス扉と外扉軽く開け、開いている状態を保ちます。

8

洗浄と消毒

コンテンツ

- ・「洗浄」上 頁 8-1
- ・「除染プロセス」上 頁 8-2
- ・「消毒と steri-run を準備する」上 頁 8-2
- ・「拭き消毒/スプレー消毒」上 頁 8-3
- ・「Steri-Run 除染モード」上 頁 8-6

洗浄

注意 利用してはいけない洗浄剤!

装置のパーツは合成物質でできています。溶剤を使うと合成物質が解ける可能性があります。強い酸やアルカリも合成物質の脆化の原因となります。合成物質製部品および表面の洗浄には、炭化水素を含んだ溶剤、10%以上アルコール含有の洗浄剤、強い酸やアルカリの入った物は使用しないでください!

水分に敏感な部品!

ディスプレイや装置背面のスイッチボックスには、洗浄剤を付けないでください。こ の部品を拭く際は、水分が入らないように気を付けてください。

外側の洗浄:

- 1。汚れの跡と堆積物を、市販の洗剤を入れたぬるま湯できちんと取り除きます。
- 2。綺麗な布で、外側を水拭きしてください。
- 3。その後乾いた布で拭いてください。

ディスプレイの洗浄:



注意水分に敏感なディスプレイ! ディスプレイは、水拭きをしたり洗浄剤を付けないでください。

・ディスプレイは 100% マイクロファイバー製の布で乾拭きしてください!

除染プロセス

Cell locker のクリーニングと消毒については、付属の Cell locker の取扱説明書を参照 してください。 装置の除染には、オペレーターが使用装置に沿った除染対策に対する衛星要綱を用意す る必要があります。 本装置には、以下の消毒方法が適切です:

拭き消毒 / スプレー消毒:

装置および補助器具の拭き消毒 / スプレー消毒は、通常の手動消毒方法です。

Steri-Run 除染モードでは、棚システムとセンサーを含む装置内空間全ての除染が自動で可能です。

消毒と ster i-run を準備する

- 1。全サンプルをサンプル室から取り出し、安全に保管します。
- 2。充分に容量のある回収容器を用意する。
- 3。ホースの先を回収容器に入れ、ホースバルブを安全に留めます。
- 4。貯水槽が空になっていく。



- 図 8-1。 給水・排水 貯水槽の排水バルブ
- 5。貯水槽の水を全て回収容器に排出し、空にする。
- 6。水の跡は布で拭きとります。
- 7。貯水槽カバーの台からエアボックス(1/図 8-2)を引き取ります。
- 8。HEPA フィルター(2/図 8-2)をエアボックスから(1/図 8-2)取り外し、エアボックスを再び取り付けます。
拭き消毒 / スプレー消毒

手動の拭き消毒 / スプレー消毒は、3 ステップあります。

- 第一消毒
- ・ 洗浄
- ・ 最終消毒

注意

・アルコール消毒剤! 10%以上アルコールを含む消毒剤は、空気に触れると発火や爆発の危険があるガ スを発生させます。

- そのような消毒剤を使用する際は、消毒プロセス中ずっと、裸火や強い熱影響を 避けてください!
- · そのような消毒剤は、よく換気がされている部屋でのみ利用すること。
- 消毒剤が効いたら、該当の装置部品を良く乾かします。
- アルコール消毒剤による発火や爆発の危険を防ぐため、安全規則に注意してください。

注意

塩化物を含む消毒剤! 塩素系消毒剤の使用禁止

注意 感電注意!

電気が通っている部分をふれると、感電し生命の危険となる場合があります。 手動の洗浄と消毒作業の前に、本装置から電力を切ってください!

- ・装置を電力スイッチで切る。
- ・ 電力プラグを抜き、誤って再接続されないようにする。
- ・装置に電力がないかを確認する。



注意 健康に危害!

装置内空間の表面は、汚染されている可能性があります。汚染された洗浄液に触れる と、伝染の可能性があります。消毒剤には、健康に危害をもたらす物質が入っている 可能性があります。

洗浄と消毒の際は、保護対策と衛生規則を守ってください!

- 保護手袋を着用。
- ・保護めがねを着用。
- ・粘膜保護のため、ロと鼻への保護も装着すること。
- 消毒剤製造者および衛生専門員からの注意に気を付けてください。

第一消毒:

1。サンプル室と取付け部品の表面は、スプレー消毒または拭き消毒を行います。



注意水分に敏感な部品!

エアダクトの裏にある CO_2 センサーと O_2/N_2 センサーには、消毒剤はかけないでください。

2。消毒剤は、製品記載のとおりに作用させてください。

取付け部品と棚システムの取り外し:

- 1。インサートプレートを取り、装置使用空間内の棚システムを全て取り外します。棚システムの取付け外しは、「棚システムを設置する」上 頁 4-8 章に記載があります。
- 2。エアボックスと HEPA フィルターを取り外します。エアボックスと HEPA フィルターの 取り付け・取り外し方法は、「HEPA フィルター 交換」上 頁 9-7 をご参照ください。



図 8-2。 HEPA フィルターとエアボックス

- 3。エアダクトの上部を(1/図 8-3)装置前方方向に引っ張り、前方の垂の鍵穴プロ フィールが装置内空間天井の留めピンをリリースしたら、下方向に浮かせて外します。
- 4。エアダクト背部の上部(2/図 8-3)を外し、装置内空間から取り出す。





- 図 8-3。 エアダクト
- 5。エアダクト背部(2/図 8-3)をから外し、取り出す。
- 6。プレフィルター(9/図 3-9)を取り出す。
 プレフィルターはオートクレーブできます。
- 7。貯水槽の蓋を取る。

装置内空間の洗浄と取外し部品:

- 1。汚れの跡と堆積物を、市販の洗剤を入れたぬるま湯できちんと取り除きます。
- 2。表面は、きれいな布で水拭きをします。
- 3。それから貯水槽から洗浄水を抜き、装置内使用空間の表面全体の水分を拭き取ります。
- 4。取外し部品を洗浄し、同様に拭いて乾かします。

最終消毒:

- 1。装置内空間の表面、棚システム、取外し部品に、改めて消毒剤をかけふき取ります。
- 2。消毒剤は、製品記載のとおりに作用させてください。
- 3。棚システムと取外し部品を再び取付けます。

Ster i-Run 除染モード

steri-runは、決まった定格温度自動の除染モードで、ヒーティング、手定格温度の維持、冷却をソフトウェアで操作します。

除染モードの全行程にかかる時間は、12時間以下です。

行程中、装置内空間では 90 分間 180 °C の乾燥した熱気が保たれ、高い除染効果があり ます。Steri-run 除染モードの効果は、独立機関によっても証明されています。ISO 11138 規格によるところの 10⁶ の菌の減少 (6 ログ減少)を達成しています。Thermo Scientific は、お問い合わせいただければこのテスト結果をお見せします。

電気機械ドアロックは steri-run 除染モード時の燃焼に対する追加の保護機能を備えて います。装置内使用空間温度が 65 °Cに達すると外扉がロックされ、危険温度中ロック され続けます。装置内使用空間温度が 65 °C以下に下がってはじめてドアロックが再び 解除されます。

行程完了後、装置を自動スタートモードで再度稼動させてください。

留意 Steri-Run 除染モードの開始妨害:

以下のエラーが検知されていると、Steri-Run 除染モードが開始できません。 ループ 温度:

- ・センサー破損、
- ・ 外扉および / または内扉(外扉が開いていると、「ドアが開いています」という エラーメッセージが表示されます)
- ・実測値が高すぎる(設定値との差異が大きすぎる)、
- ・実測値が低すぎる(設定値との差異が大きすぎる)、
- ・実測値 妥当ではない、
- 接続障害。
- ・水が検出

過温度保護機能:

装置の過温度保護機能がオンになると、Ster i-Run 除染モードはエラーが解消した後に 開始することができます。

Ster i-Run 除染モード中にガス供給がない、ガスモニタリングシステム(オプション) 付き装備の場合:

Steri-Run 除染モード中にエラー 「ガスの不足」が表示されると、聴覚シグナル (ブ ザー音)が鳴ります。このシグナルは、ディスプレイに触れると鳴りやみます。この場 合、Steri-Run 除染モードは停止しません。アラームリレーは、ガスモニターが満タン のボトルを認識するまで、作動したままとなります。



注意 Cell locker の使用耐熱温度! 「除染実行」の前に、Cell locker を取り出す必要があります。使用耐熱温度は最高 121°C (250°F)です。 Ster i-Run 除染モードのプロセス:

- 1。除染モードを開始する前に、内側容器のシリコンストッパーを取り外し、通路開口 部の外側に差し込んでください。
- 2。棚システムの部品は、洗浄の後に装置内空間に再び取り付けます。
- 3。装置を電力スイッチでオンにする。
- 4。除染モードをオンにし、開始します。
- 5。Steri-run 運転完了後、装置のスイッチを切ってください。
- 6。エアボックス (1/図 8-2) を取り外し、HEPA フィルター を再度取り付けます (2/図 8-2)。
- 7。必要に応じて、自動スタートで運転を再開してください。



ガラス扉の取手とガラスは、外扉インナーパネルおよび表面棚システムと装置内空間 は、Steri-Run 除染モード中大変熱くなります。

モードの運転中もしくは中断後すぐはこれらの面には保護手袋を装着してから触れて ください!

ドアロック・キット付のインキュベーターは、Steri-run 除染モードによって内部表面が高温の間、装置内へのアクセスを防ぎます。

注意 サンプルの破損!



装置内空間は、Steri-Run 除染モード 中は 180 °C に熱されています。以下を確認し てください:

- · 全サンプルが装置内空間から取り出されたか、
- ・ 補助器具は全て装置内空間から取り出されたか。

Ster i-Run 除染モード運転プロセス:

Steri-Run 除染モードの残り時間は、開始もしくは現時点から冷却プロセスが終わるまでの時間です。表示されている残り時間は計測値ではなく、目安の値です。 モードは 3 つの プロセスから成っています:

- 1。ヒーティングプロセス、
- 2。除染プロセス、
- 3。冷却。

ヒーティングプロセス:約2h 装置内空間は180°Cに熱されます。 電磁性ドアロック・キット装備のインキュベーターは、装置内部温度が65°Cを超過す るとドアをロックします。

除染プロセス: 約 1.5 h 除染温度環境を作り出した後、約 90 分の除染プロセスが開始します。その際温度は 180 °C に保たれます。 冷却プロセス:約8h 本装置は元々設定されている温度設定値まで冷却します。 電磁性ドアロック・キット装備のインキュベーターは、装置内部温度が65°C以下に下 がるドアロックを解除します。



図 8-4。 除染モードのプロセス

Steri-Run をオンにする

Ster i-Run は、自動の除染モードで、装置内空間の消毒を行います。

1。Steri-Run キーを押します。

留意

シリコンの磨耗を防ぐため、Steri-run 除染を始める前に、内側容器から必ずシリコンストッパーを取り外し、通路開口部の外側に差し込んでください。

・メニュー Steri-Run 操作方法が表示されます。



図 8-5。 メニュー Steri-Run 操作方法

- 2。メニュー Steri-Run 操作方法を終了し、Steri-Run を中断する:
- · 終了キー を押します。
- ・メインメニューに戻ります。

- 3。Steri-Run をオンにする:
- · START ボタンで起動します。
- · Steri-Run プロセスのダイアログボックスが表示されます。
- 4。時間ブザーが鳴るまで 30 秒間、装置内使用空間を換気し、装置のドアを両方とも開 けます。
- 5。全サンプルを装置内空間から取り出します。
- 6。貯水槽の排水をし、水滴を拭き取ります。
- 7。時間シグナルが鳴ったら、装置ドアを両方とも閉めます。
- Steri-Run を開始する。
- Steri-Run 除染モードのプロセスが開始します。装置が温まり、65°C で電気機械
 ドアロックが作動します。
- Ster i-Run 除染モード運転中は、ディスプレイに現在のステータスが表示され、以下の情報が示されます:
 - ・温度,
 - · 開始時間、
 - ・プロセス、
 - ・残り時間。

Start	steri-run - ステータス 温度: 57 ℃ 開始時間: 13:42 プロセス: 停止 残り時間: 07:01
	Stop

図 8-6。 Steri-Run をオンにする

Steri-Run を中断する

Steri-Run 除染モードは、いつでも中断させることができます。

- 1。Steri-Run を中断する:
- · Stop キーを押します。
 - Stop ボタンを押すと、安全データ確認としてダイアログボックス steri-run Stop が呼び出されます。ここで完全に停止させてしまうことも、再開すること もできます。
- 2。Steri-Run を終了:
- ·終了キーを押します。

- エラーが表示されます。
- · エラー表示を確定すると、メインメニューに戻ります。
- 3。Steri-Run を続行する:
- · **戻る**キーを押します。
- · ステータス表示へ戻り、除染モード が続行されます。
- 4。Steri-Run をステータス表示画面から中断させる:
- Stop キーを押します。
- 安全データ確認として、ダイアログボックス ster i-run Stop が呼び出されます。続きはステップ2へ。

Steri-Run のエラー中断

除染モード運転中にエラーが発生すると、エラーが表示され以下の行動に誘導されます:

- ・ 除染モードが自動的に冷却プロセスに変わる。
- ・聴覚シグナル(ブザー音)が鳴る。
- 1。聴覚シグナルを確定する:
- ディスプレイに触れます。
- ・聴覚シグナルが鳴りやみます。終了キーが表示されます。除染モードを手動でキャンセルしない場合は、設定温度まで冷却され、装置内が乾燥されます。
- 2。Steri-Run を中断する:
- · 終了キー を押します。
- エラーが表示されます。
- エラー表示を確定すると、メインメニューに戻ります。

S S	teri-run 終了> 消 戻る> 継	- 中 i毒プ 続す	断 ロセスの停」 る	Ł	
	₽.Z		(1) (1) (1) (1) (1) (1) (1) (1) (1) (1)	<u> </u>	

図 8-7。 Steri-Run を停止、中断する

Steri-Run 終える

これらの3段階が完全に完了すると、自動的にダイアログボックス steri-run 完了 (図 8-8)が表示されます。除染モードは手動で終える必要があります。

steri-run - 終了 1.ドアを開く 2.水を補充する 3.HEPAフィルターを取付ける 4.ドアを閉める 5.必要に広じて自動スタートモードを実行
5.必要に応じて自動スタートモートを実行

図 8-8。 Steri-Run を終了

- ・ Steri-Run を終了:
 - ·終了 キー を押します。
 - ・メインメニューに戻ります。

留意 Steri-run 除染モード中にドアを開ける:

除染モード中にドアを開け閉めすると、そのプロセスを最初からやり直しし、モード がエラーなく続行するようになっています。

留意

燃焼の危険性!緊急の際を除きドアを開けないでください。

Steri-run 除染モード中断後にドアを開ける

電気機械ドアロック・キット付の装置のドアは、Steri-run 除染モード中断直後はユー ザーあるいはエラーによって開けることができません。

負傷しない温度の 65 °C に下がる前に外側ドアのロックを解除するためには、緊急用始 動レバーを引く必要があります。(図 8-9 の Pos. 3):

・ 緊急用始動レバー(3)の位置をインキュベーターの底部に合わせます。レバーは装置左側面、電力スイッチ(4)から約5 cm 横/後にあります(図 8-9 下参照)。



図 8-9。 インキュベーター下部のドアロックレバーと緊急用ロック解除

・ 底面から手を伸ばして緊急用始動レバー(3)を引きます。

ドアロックが解除されました。外側ドアを開けることができます。



注意 表面熱い!

ガラス扉の取手とガラスは、外扉インナーパネルおよび表面棚システムと装置内空間は、Steri-Run 除染モード中大変熱くなります。

モードの運転中もしくは中断後すぐはこれらの面には保護手袋を装着してから触れて ください!

9

メンテナンス

コンテンツ

- ・「検査とコントロール」上 頁 9-1
- ・「メンテナンスの間隔」上 頁 9-2
- ・「温度キャリブレーションの準備」上 頁 9-2
- ・「温度の整合を行う」上 頁 9-3
- ・「CO2 キャリブレーションを準備する」上 頁 9-5
- 「CO2 キャリブレーションを行う」上頁 9-6
- ・「HEPA フィルター 交換」上 頁 9-7
- ・「ガスダクトフィルター交換」上 頁 9-9
- 「装置安全装置の交換」上 頁 9-10
- ・「ドアパッキンの交換」上 頁 9-10

検査とコントロール

装置の機能と安全性を保つため、以下の機能と装置部品の検査は、それぞれ一定の期間で もって行ってください。

Cell locker のメンテナンスについては、付属の Cell locker 取扱説明書を参照してください。

- 毎日のコントロール:
 - · CO₂ 供給設備のガス量
 - 0₂/N₂ 供給設備のガス量

年次検査:

- ・ ガラス扉パッキンの密閉具合
- ・ 圧力調整口の流れ具合
- ・コントロールパネルと装置調整の操作テスト
- ・ 電気安全検査は、国内の規約に則ります(例 BGV 3)

留意 機能の確認:

検査のために安全装置が取り外されたり機能しない状態になる場合、安全装置が再び 取り付けられ問題なく機能するか確認がされたて初めて、本装置を再び使用すること ができます。

メンテナンスの間隔

装置使用期間中は、以下のメンテナンスを実施してください:

四半期毎のメンテナンス:

・ 自動スタートモードと Ster i-Run 除染モードを運転する。

・ 温度と CO₂/O₂ 比較計測を実施する。

6分割ガスシェード:

・カバーガスケットは6か月ごとに確認します。必要であれば、交換します。

・メンブレンフィルターについては、付属の取扱説明書を参照してください。

年次メンテナンス:

・ガスダクトフィルターを交換する。

・ テクニカルサービスのサービス検査を実施する

留意メンテナンス契約:

Thermo Scientificは、装置に合ったメンテナンスの契約をご提供いたします。この契約には、必要な検査と手入れが含まれています。

温度キャリブレーションの準備

装置内部温度センサーが正確な計測値を示すために、四半期に一度温度の比較計測を実施してください。その際に計測に大きな差異が確認された場合、温度整合を行ってください。その際、装置の温度調整は比較計測の計測値に設定されます。 比較計測には、正確性 < ± 0.1 °Cに校正された計測器を使用してください。 計測中の温度変化を最小限に抑えるため、計測器は等温容器(例えばグリセリンの入った容器など)に入れて、装置内空間に置いてください。比較計測は、装置内空間中央部で行います。

留意 等温容器: 等温容器として、水で満たされた容器は使わないでください。水の蒸発により、低す ぎる温度が算出されてしまいます。

装置内空間の温度が高すぎる: 整合後に装置内空間温度が上がってしまうことがあるので、扉を約30秒間開けて解 消してください。



図 9-1。 温度キャリブレーションの準備

比較計測を実行する:

- 1。装置を電力スイッチでオンにする。
- 2。 温度設定値を設定し、本装置がウォームアップされるのを待つ。これには、数時間を 要する場合があります。
- 計測器(2)を装置内空間中央のインサートプレートの中心に置く。
 他の方法としては、温度センサーと同じポジションでの配置も可能です。管はガラス 扉の計測口(3)または装置背面のカバーロゼット(1)に移されます。
- 4。扉を締める。
- 5。計測器の温度が安定するまで待つ。
- 6。温度の整合を行う

温度の整合を行う

計測例:

- ・ 温度の設定値 t: 37 ° C
 計測した比較温度: 36.4 ° C
- 1。温度表示 キー を押します。 ・温度メニュー (図 9-2)が表示されます。



図 9-2。 温度表示と温度選択メニュー

- 2。温度メニューを終了する:
- · 終了キー を押します。
- 3。サブメニュー 校正値を呼び出す:
 校正キー を押します。



図 9-3。 温度の整合を行う

4。計測した値(ターゲット値)を入力する: 時間値は、キーを押すことで段階的に高くしたり低くしたりすることができます。-キーまたは+キー長押しをすることで、数値は高速で変化していき、約3秒間長押し を続けると、さらにその速度がアップします。

ターゲット値を上げる: ・+キーを押します。

設定値 を下げる:

- ・ -キーを押します。
- 5。ターゲット値を 適用し保存する:
- · Enter **キー** を押します。
- ・ **保存**キーを押します。

・メインメニューに戻ります。装置内空間にて計測された現在の実測値が、温度表示
 部分に表示されます。

留意 装置内空間の温度が高すぎる:

整合後に装置内空間温度が上がってしまうことがあるので、扉を約 30 秒間開けて解 消してください。

値を戻す:

30 秒間値の変更が行われない場合、変更メニューは自動的に終了し、最後に保存された 値のままとなります。

CO₂ キャリブレーションを準備する

装置内部 CO₂ センサー計測値の的確さ保つため、四半期に一度 CO₂ 比較計測を行ってくだ さい。

その際に計測に大きな差異が確認された場合は、 CO_2 キャリブレーションを行ってください。 その際、装置の CO_2 調整は比較計測の計測値に設定されます。比較計測には、正確性 $< \pm 0,3\%$ CO_2 に校正された計測器を使用してください。

適切な計測器具:

- 持ち運び可能な IR ポータブル計測器具。
 - 計測サンプルは、閉めることが可能なガラス扉の計測口から取ります。比較計測は、 ウォームアップした装置にて行ってください。

比較計測を実行する:

- 1。装置を電力スイッチでオンにする。
- 2。CO2の設定値を設定し、装置がウォームアップされ湿度環境を構築するのを待ちます。
- 3。IR ポータブル計測器具の測定プローブは、計測口から装置空間に入れます。計測器 具が安定した CO₂ 値を出すまで待ちます。
- 4。オプションのガスシェード搭載の装置では、計測口は以下のところにあります:
- STERI-CYCLE i160 LK では中央ガスシェード、
- ・ *STERI-CYCLE i250 LK* では左中央ガスシェード。
- 5。測定プローブを抜き、計測口を閉めて扉も閉じます。
- 6。CO₂ 調整を校正する。

留意 IR 測定セル: IR 測定セル付き装置の場合、CO₂ キャリブレーションは CO₂ 濃度が CO₂ 4.0%以上に 設定されている場合に、行うことができます。 校正は、作業プロセスで予定している CO₂ 設定値(将来の作業プロセス値)にて実施 してください。



図 9-4。 ガスシェードの計測口

CO₂ キャリブレーションを行う

計測例: ・ CO₂ 設定値: 5 % サンプル計測: 5.6 %

1。CO₂ 表示キーを押します。 · CO₂ メニューが表示されます。



- 2。CO₂メニューを終了する:
- · 終了キー を押します。
- 3。サブメニュー 校正値を呼び出す: ・ 校正キー を押します。



- 図 9-5。 CO₂ キャリブレーションを行う
- 4。計測した値(ターゲット値)を入力する: 時間値は、キーを押すことで段階的に高くしたり低くしたりすることができます。-キーまたは+キー長押しをすることで、数値は高速で変化していき、約3秒間長押し を続けると、さらにその速度がアップします。

ターゲット値を上げる:

・ + キーを押します。

設定値 を下げる:

- ・ ーキーを押します。
- 5。ターゲット値を 適用し保存する:
- · Enter **キー** を押します。
- · 保存キーを押します。
- ・メインメニューに戻ります。装置内空間にて計測された実測値が、CO₂表示部分に表示されます。

留意 CO₂ 含有量が高すぎる: 整合後に CO₂ 含有量が上がってしまうことがあるので、扉を約 30 秒間開けて解消して ください。

値を戻す:

30 秒間値の変更が行われない場合、変更メニューは自動的に終了し、最後に保存された値のままとなります。

HEPA フィルター 交換

HEPA フィルターは、装置空間背面前の底にあるプラスチックカバー(エアボックス)の下にあります。

HEPA フィルターの交換作業工程:

- 1。装置をオフにし、ガス供給を停止して、装置内空間を換気する。
- 2。中央と下のインサートプレートを装置内空間から取り出す。



3。貯水槽カバーの前面を上にあげます(1/図 9-6)。

- 図 9-6。 エアボックスを取り外す
- 4。貯水槽カバーを前に引っ張ります(2/図 9-6)。
- 5。HEPA フィルター付エアボックスを取り外します。
- 6。エアボックス(1/図 9-7)を回し、エアボックスの左側の垂 を HEPA フィルターのかみ合い(6/図 9-7)から外します。
- 7。エアボックス(1/ 図 9-7)右側の垂(3/ 図 9-7)を HEPA フィルターの該当の溝(4/ 図 9-7)から引き抜きます。









図 9-7。 HEPA フィルターの取付け

- 8。新しい HEPA フィルターをエアボックス(1/図 9-7)に取り付け、留めます。
- 9。エアボックスを貯水槽カバーの台の上に設置します。
- 10。装置が HEPA フィルターなしで運転されたことがある場合、説明書通りに 「HEPA フィ ルターをオン / オフにする:」上 頁 6-39 ユーザーコンフィギュレーションにて HEPA フィルターをオンにしてください。
- 11。必要に応じて、説明書に沿って「リマインダーを設定する」上 頁 6-26 ユーザーコンフィギュレーションにて HEPA フィルター交換のリマインダーを設定します。HEPA フィルター交換のリマインダーは、1~12か月の中で設定ができます。設定値は、発送時の初期設定である6か月を奨励しています。

ガスダクトフィルター交換

ガスダクトフィルター(CO₂/O₂/N₂供給)は、プラスチックねじが付いており、スイッチ ボックスわきのネジのレセプタクルにしっかりと留めます。

ガス供給のガスダクトフィルター交換作業工程:

- 1。ガス供給が閉じていることを確認する。
- 2。ホースクランプ(4/図 9-8)を緩めます。

3。ガスホース(5/図 9-8)をガスダクトフィルターの接続ノズルから引き離します。

全ガスダクトフィルターの作業方法:

- 4。セーフティプレート(1/図 9-8)のネジを外します。
- 5。ガスダクトフィルター(2/図 9-8)を、ネジのレセプタクル(3/図 9-8)から取り外 します。
- 6。ガスダクトフィルターを回しはめる際は、プラスチックねじがクロススレッドしてい ないよう気を付けます。ガスダクトフィルターを注意深く回しはめる。
- 7。セーフティプレートをねじで固定する。

ガス供給のガスダクトフィルター交換作業工程:

8。ガスホースをフィルターの接続ノズルに乗せ、ホースクランプで固定します。ガス ホースが接続ノズルにきちっと装着されているかを確認してください。



図 9-8。 ガスダクトフィルター 取付け

装置安全装置の交換

安全装置の交換は、装置利用者が行うことができません。技術的エラーによって装置が 停止してしまった場合、テクニカルサービスを呼んでください。

ドアパッキンの交換

留意

扉のパッキンの交換は、テクニカルサービス員か資格のある専門員に頼むことをお勧めいたします。



コンテンツ

·「使用素材一覧:」上 頁 10-1

注意 汚染の危険性があります!



本装置は、感染物質の処理に使われていた可能性があります。それゆえ、本装置および装置の部品は、汚染されている可能性があります。装置の構成部品は、廃棄処分前 に必ず除染すること!

- ・装置の構成部品はしっかりと洗浄し、使用用途によっては更に殺菌や滅菌を行っ てください。
- ・廃棄処分品には、除染処置実行に関する明確な記載をした非危険物証明書を添付 すること。

装置の部品は全て、決められた廃棄処分方で除染した上で、引き渡しが可能です。 HEPA フィルターは、その国内で有効な特別廃棄物の規定に沿って廃棄すること。

留意 リサイクルサービス: Thermo Fisher Scientific社は、環境に配慮した使用済み装置のリサイクルサービス をご提供しています。

使用素材一覧:

部品	素材
断熱部品	グラスウール、片面被覆グラスフリース付 グラスウール
電子基板	様々な合成物質でカバーしてある電子部 品、エポキシ樹脂の基板上に装備
合成物質の部品全般	ABS 樹脂および PPS 樹脂 GF40、素材ラベ ルにご注意下さい
外装	亜鉛メッキをしラッカー塗料した薄鋼板、 ステンレス鋼 1.4016
装置背面	亜鉛メッキ加工薄鋼板

部品	素材
外扉	亜鉛メッキをしラッカー塗料した薄鋼板、 ステンレス鋼 1.4016
ドアインナーパネル	ステンレス薄鋼板 1.4301
操作および表示ディスプレイフィルム	ポリエチレン
ヒーター	シリコン被覆の耐熱導体
内側容器、内蔵品、インサートプレート	ステンレス鋼 1.4301、 銅
圧カ調整ロへの使用	ステンレス鋼 1.4301(ソケット)、 1.4404(焼結フィルター)
ガラス板	ケイ酸ガラス
センサーブロック (WLD)	ステンレス鋼 1.4301
ケーブル	合成物質およびシリコン被覆の銅線
エラストマ、全般	シリコン
フィルター	HEPA フィルター マイクロファインガラス, ABS 樹脂・シリコン製 Cell locker 用メン ブレンフィルター、 ガスフィルター ポリプロピレン容器と GF/PTFE 膜、 プレフィルター 金属布 ステンレス鋼 1.4401
包装	段ボール、ポリエチレンシート、発泡スチ ロール成形品、ポリプロピレン
ドアパッキン用マグネット	永久磁石
Cell Locker	ポリカーボネート製 Makrolon 2528

仕様

コンテンツ

- ・「STERI-CYCLE i160 LK」上 頁 11-2
- ・「STERI-CYCLE i250 LK」上 頁 11-6

11

STERI-CYCLE i160 LK

名称	ユニット	値
機械		
外寸(幅 x 高さ x 奥行)	mm	637 x 905 x 790
内寸(幅 x 高さ x 奥行)	mm	470 x 607 x 576
内部容量、 うち使用容量		約 165 約 100
インサートプレート(幅 x 奥行) 数、納品量 数、最大 分布荷重、最大	mm 個 個 kg	423 x 465 3 11 10 / インサートプレート
装置合計荷重、最大	kg	30
重量	kg	83
重量、付属品含まず	kg	78. 4
<u>熱</u>		
DIN 規格の熱感知安全装置 12880:2007-05		クラス 3.1 (超過温度の場合の監視機能 付き温度監視機 (TWW))
周囲温度域	°C	+1834
積重ねた装置の周辺温度	°C	+1828
温度調整域、インキュベーション	°C	RT + 355
時間的な温度差異(DIN 12880, パート 2)37 °C時	°C	± 0.1
空間的な温度差異(DIN 12880、パート 2) 37 °C時 ^{*1})	°C	<= ± 0.3
自動スタートモードにかかる時間: 37 °C時 周辺温度 20 °C	h	5 10
周囲への熱出力: 37 °C時 Steri-Run 除染中	kWh/h kWh/h	0. 06 0. 59
湿度		
水質		電気抵抗: 50 kOhmcm ~ 1 MOhmcm 伝導率: 1 ~ 20 μS/cm
補填量 [:] インキュベーション運転	I	最大 3 / 分 0.5
37 °C時の持続湿度(高湿度モード) 37 °C時の持続湿度(低湿度モード)	%rH %rH	約 93 約 90

名称	ユニット	値
その他		
音圧レベル (DIN 45 635、パート 1)	dB(A)	< 50
周囲の相対湿度	%rH	最大 80
設置場所の高さ	m NN	最大 2000

*¹⁾ 装置通常運転における DIN 12880 に則った数値 校正仕様に記載のある値を採用すること

STERI-CYCLE i160 LK

名称	ユニット	值
C0 ₂ ガス技術		
ガス純度	%	最低 99.5 または医療品質
見本	bar	最低 0.8 ~最大 1 (11.6 - 14.5 psi)
測定・調整範囲	Vol %	0 20
時間的な調整差異	Vol %	± 0.1
CO ₂ 測定セル		
37 °C で CO ₂ が 5% 時の正確性	% CO ₂	± 0.3
0 ₂ ガス技術		
ガス純度	%	最低 99.5 または医療品質
見本	bar	最低 0.8 ~最大 1 (11.6 - 14.5 psi)
測定・調整範囲	Vol %	1 21 または 590
時間的な調整差異	Vol %	± 0.2
02 測定セル		
37 °Cで0 ₂ が21%時の正確性	%0 ₂	± 0.5 (オプション: 121% 0 ₂) ± 2.0 (オプション: 590% 0 ₂)
電子的		
定格電圧	V	$\begin{array}{cccccccccccccccccccccccccccccccccccc$
定格周波数	Hz	50/60
保護方 (IEC 60529)		IP 20
保護クラス		I
過電圧カテゴリ(EN 61010)		II
汚染度(EN 61010)		2

名称	ユニット	值
定格電流	A	230 V: 除染: 4.6 インキュベーション: 2.4 220 V: 除染: 4.4 インキュベーション: 2.3 120 V: 除染: 8.3; インキュベーション: 4.6 100 V: 除染: 7.2; インキュベーション: 3.9
導線保護スイッチ		16 A
定格消費電力	kW	230 V: 除染: 1.10 インキュベーション: 0.56 220 V: 除染: 0.97 インキュベーション: 0.51 120 V: 除染: 1.01 インキュベーション: 0.55 100 V: 除染: 0.72 インキュベーション: 0.39
EMV クラス		В

STERI-CYCLE i250 LK

名称	ユニット	值
機械		
外寸(幅 x 高さ x 奥行)	mm	780 x 970 x 945
内寸(幅 x 高さ x 奥行)	mm	607 x 670 x 629
内部容量、 うち使用容量	Ι	約 255 約 162
インサートプレート(幅 x 奥行) 数、納品量 数、最大 分布荷重、最大	mm 個 個 kg	560 x 500 3 12 銅製シェルフあたり 10 ステンレススチール製シェ ルフあたり 14
装置合計荷重、最大	kg	30 銅製シェルフ/ 42 ステンレススチール製 シェルフ
重量、付属品含まず	kg	97. 5
熱		
DIN 規格の熱感知安全装置 12880:2007-05		クラス 3.1 (超過温度の場合の監視機能 付き温度監視機 (TWW))
周囲温度域	°C	+1834
積重ねた装置の周辺温度	°C	+1828
温度調整域	°C	RT + 355
時間的な温度差異(DIN 12880、 パート 2)	°C	± 0.1
空間的な温度差異(DIN 12880、パート 2) 37 °C時 ^{*1)}	°C	± 0.3
自動スタートモードにかかる時間: 37 °C時 周辺温度 20 °C	h	5 10
周囲への熱出力: 37 °C時 Steri-Run 除染中	kWh/h kWh/h	0. 07 0. 75
湿度		
水質		電気抵抗: 50 kOhmcm ~ 1 MOhmcm 伝導率:1 ~ 20 μS/cm
補填量: インキュベーション運転		最大 3 / 分 0.5

名称	ユニット	値
37 °C時の持続湿度(高湿度モード)	%rH	約 93
37 °C時の持続湿度(低湿度モード)	%rH	約 90

その他		
音圧レベル (DIN 45 635、パート 1)	dB(A)	< 50
周囲の相対湿度	%rH	最大 80
設置場所の高さ	m NN	最大 2000

*1) 装置通常運転における DIN 12880 に則った数値 校正仕様に記載のある値を採用すること

STERI-CYCLE i250 LK

名称	ユニット	值
C0 ₂ ガス技術		
ガス純度	%	最低 99.5 または医療品質
見本	bar	最低 0.8 ~最大 1 (11.6 - 14.5 psi)
測定・調整範囲	Vol %	0 20
時間的な調整差異	Vol %	± 0.1
CO ₂ 測定セル		
37 °C で CO ₂ が 5% 時の正確性	% CO ₂	± 0.3
0 ₂ ガス技術		
ガス純度	%	最低 99.5 または医療品質
見本	bar	最低 0.8 ~最大 1 (11.6 - 14.5 psi)
測定・調整範囲	VoI %	1 21 または 590
時間的な調整差異	VoI %	± 0.2
02 測定セル		
37 °Cで0 _{2が21%} 時の正確性	%0 ₂	± 0.5 (オプション: 121% 0 ₂) ± 2.0 (オプション: 590% 0 ₂)
電子的		
定格電圧	V	$\begin{array}{cccccccccccccccccccccccccccccccccccc$
定格周波数	Hz	50/60
保護方 (IEC 60529)		IP 20
保護クラス		Ι
過電圧カテゴリ(EN 61010)		II
汚染度(EN 61010)		2

名称	ユニット	值
定格電流	A	230 V: 除染: 5.5 インキュベーション: 3.3 220 V: 除染: 5.3 インキュベーション: 3.2 120 V: 除染: 10.4
		インキュベーション: 6.3 100 V: 除染: 8.9 インキュベーション: 5.3
導線保護スイッチ		16 A
定格消費電力	kW	230 V: 除染: 1.26 インキュベーション: 0.76 220 V: 除染: 1.16 インキュベーション: 0.69 120 V: 除染: 1.25 インキュベーション: 0.75 100 V: 除染: 0.89 インキュベーション: 0.53
EMV クラス		В

11 仕様 STERI-CYCLE i250 LK

添付物 データ通信

コンテンツ

- ・「データコミュニケーションのコマンドシーケンス構造」上 頁 12-4
- ・「一般パラメータの一覧 (アドレス Oxxx)」上 頁 12-5
- ・「一覧インキュベータ・パラメータ (アドレス n 2xxx)」上 頁 12-5
- ・「構造 エラーメモリ」上 頁 12-8
- ・「データログ構造」上 頁 12-11
- 「コード例 データログ」上 頁 12-15
- ・「プログラム STERI-CYCLE i160 LK / STERI-CYCLE i250 LK」上 頁 12-21

USB インターフェース

装置には USB インターフェースが一つ搭載されています。USB インターフェースは、USB 1.1 / USB 2.0 / USB 3.0 (フルスピード)のスタンダードに適合します)。USB インター フェースは、仮の Com ポートとして動きます。それゆえ、インターフェースの伝達速度 を定義されたボーレート (9,600、19,200、38,400、57,600 ボー)内に変更することが できます。データ交換は定められたコマンドシーケンス構造を通じて行われます。コマ ンドシーケンスは RS 232 インターフェースの構造パターンに則ります。

留意 仮 Com ポート付の USB 接続の設定:

パソコンとインキュベータ間のデータ交換に USB インターフェースを使用する場合 は、納品時同封の仮の Com ポート(USB シリアルポート)としての USB 接続のドライ バを使用し、設定を行います。 装置管理・接続の Windows ダイアログボックスにて、Com ポートの割り当てがわかり ます。例: USB シリアルポート(COM5)にはプログラム *STERI-CYCLE i160 LK / STERI-CYCLE i250 LK* が接続インターフェースとして定められます(参照 「プログラム STERI-CYCLE i160 LK / STERI-CYCLE i250 LK」上 頁 12-21)。

🗏 Device Manager	
<u>Eile A</u> ction ⊻iew <u>H</u> elp	
🖶 🕮 Network adapters	*
🖶 🖉 Ports (COM & LPT)	_
- 🖉 Communications Port (COM1)	
- 🦉 Communications Port (COM2)	
- 🖉 ECP Printer Port (LPT1)	
- 🖉 USB Serial Port (COM3)	
🕀 🚓 Processors	=
🗄 🧐 Sound, video and game controllers	
🖶 😼 System devices	
🗄 🗳 Universal Corial Due controllere	<u> </u>

ドライバは、以下の OS に対応しています: WIN 7, WIN 8, WIN 2000, WIN XP, WIN VISTA。

ドライバ USB インターフェースをインストールする

USB ケーブルをスイッチボックス **STERI-CYCLE i160 LK** / **STERI-CYCLE i250 LK** の所にある USB イ ンターフェース用(オプション)に接続し、パソコンと繋げます。 Windows のハードウェア探知が USB 接続を認識したら、「新しいハードウェアの検索ウィ ザード」のダイアログボックスが開きます。

1。「ソフトウェアの検出はしない」というオプションを選択。



2。「ファイル名を特定しソフトウェアをインストールする」を選択。

Assistent für das Suchen neuer Hardware	
	Mit diesem Assistenten können Sie Software für die folgende Hardwarekomponente installieren: EVAL232 Board USB <> Serial Falls die Hardwarekomponente mit einer CD oder Diskette geliefert wurde, legen Sie diese jetzt ein.
	Wie möchten Sie vorgehen?
	 ○ Software <u>a</u>utomatisch installieren (empfohlen) ○ Software von einer Liste oder bestimmten Quelle Installieren für fordreschöttene Benutzen)
	Klicken Sie auf "Weiter", um den Vorgang fortzusetzen.
	< Zurück Weiter > Abbrechen

3。CD-ROM を指定。



4。CD-ROM内でドライバを選択します。

Browse For Folder	?×
Select the folder that contains drivers for your hardware.	
🗉 🛗 My Documents	~
🗉 😼 PC15	-
Image: Bigging (A:) Image: Bigging (A:)	
🗉 🥯 Festplatte_PC15 (C:)	
🛛 🥝 HERAcell_v01 (D:)	-
🖃 🗁 _Driver	-
🗀 amd64	
i386	
🗀 _Programs	
DE DE	
ES ES	
E FR	-
🗋 GB	*
Ta ulau anu subfaldera, slista alus sina sharus	
to view any subrouers, click a plus sign above.	
	ancel
	.:

5。インストールウィザードが以下のドライバのインストールを行います: EVAL22 Board USB インストールが完了したら、ウィザードを閉じてください。

インターフェースの伝達速度は、定義されたボーレート(9,600、19,200、38,400、 57,600 ボー)内においてタッチディスプレイ **STERI-CYCLE i160 LK** / **STERI-CYCLE i250 LK** にて設定ができます(参照「設定 / セットアップ」上 頁 6-17)。

データコミュニケーションのコマンドシーケンス構造

パソコンとインキュベーター STERI-CYCLE i160 LK / STERI-CYCLE i250 LK 間でのデータのやり取 りにおいて送受信したデータ記号は、全て ASCII コードで、通常のターミナルにおいて 表示可能です。 これにより、装置使用の開始、コミュニケーション管理とプログラミングを、簡単に行 うことができます。

記録の説明

記号のコーデック: ASCII 記号、大文字は使用不可能。

パラメータを読み取る:

?:aaaa:bb::cc <cr></cr>
?:aaaa:bb:XXXX:cc <cr></cr>
!:aaaa:bb:XXXXX:cc <cr></cr>
aaaa = パラメータのアドレス
bb =該当テレグラムにおける使用データの数(00 - ff)
cc = チェック数: CRC8-CCITT: x8 + x2 + x1 + 1 = 0x07
cc なしおよび〈CR〉
XXXX = bb バイトの使用データ

返答要素の説明:

aaaa	パラメータのアドレス
bb	該当テレグラムにおける使用データの数(00 - ff)
cc	チェック数: 全バイトの投入 XOR
	チェック数なし <cr></cr>

ソフトウェアバージョンデータ確認の例(50111927): データ確認: ?:0001:00::cc<CR>

返答: !:0001:08:50111927:cc<CR>

パラメータを書く:

コマンド: !:aaaa:bb:XXXXX:cc<CR>

- 返答: !:aaaa:bb::cc<CR>
- 説明: aaaa = パラメータのアドレス bb = 該当テレグラムにおける使用データの数(00 - ff) cc = チェック数: CRC8-CCITT: x8 + x2 + x1 + 1 = 0x07 cc なしおよび <CR> XXXX = bb バイトの使用データ

エラー表示のある返答:

返答: !:aaaa:bb:XX:cc<CR>
返答要素の説明:
 aaaa パラメータのアドレス,
 bb 使用データの数(常に 02)
 cc チェック数: CRC8-CCITT: x8 + x2 + x1 + 1 = 0x07
 cc なしおよび <CR>
 XX = 2 バイト エラー表示(下記の表参照)

不明コマンドの例:

問合せ:	?:0005:00::cc <cr></cr>

返答!:0005:02:?1:cc<CR>

エラー表示における両バイトの意味:

エラー表示	説明
?0	テレグラム構造またはチェック数におけるエラー
?1	不明コマンドまたは不明パラメータ
?2	内部保存エラー
?3	データエラー(値は限界値の範囲内)

ー般パラメータの一覧 (アドレス Oxxx)

一般パラメータは、日付、時刻、メインプレートのバージョン番号などのシステム値です。

一般パラメータを読み取る

アドレス	説明	注
0001	バージョン番号 メインボード	8 桁
0010	アウトプット 日付と時刻 [時:分:秒]; [日:月:年]	17 バイト / 十進値 フォーマット xx:xx:xx;xx:xx:xx
0011	日付 [日︰月︰年]	8 バイト / 十進値 フォーマット xx:xx:xx
0012	時刻 [時:分:秒];	8 バイト / 十進値 フォーマット xx:xx:xx

ー覧インキュベータ・パラメータ (アドレス n 2xxx)

インキュベータ・パラメータは、以下の通りに区分されています:

・3つのループ温度のパラメータ (Basic)、 CO_2 、 O_2 、

・オペレーティング機能とデータ記録のパラメータ (内部機能)

パラメータ(Basic)を読み取る

アドレス	説明	注
2000	装置ステータス ^{*1)} (エラーおよび) ループのステータス 温度, C0 ₂ 、0 ₂ 、rH、参考温度。	33 バイト / 十六進値 フォーマット xxxxxxx;xxx; ;xxxx;xxxx;xxxx
2010	設定、実測、参考温度 ^{*2})	23 バイト / 十進値 フォーマット +xxx. xx;+xxx. xx;+xxx. xx
2020	設定および実測 CO ₂ 含有量 ^{*2)}	15 バイト / 十進値 フォーマット +xxx.xx;+xxx.xx
2030	設定・実測 0 ₂ 含有量 ^{*2)}	15 バイト / 十進値 フォーマット +xxx.xx;+xxx.xx
204a	実測水位(100% または 0%)	7 バイト / 十進値 フォーマット +xxx.xx
204b	ディスプレイ 低湿度(1 オン, 0 オフ)	2 バイト / 十六進 フォーマット xx

*¹⁾ 例 装置ステータスおよび(エラーの)ループステータス (詳細はエラー表示の表を参照)

*2) 全値は、小数点第二位までの数字です。

パラメータ(内部機能)を読み取る

アドレス	説明	注
2100	ステータス 経過 ^{*1)} と 残り時間 [時感 [:] 分] 消毒、および前の開始日付 と時刻	25 バイト / 十進値 フォーマット xx;+xxx:xx;xx.xx.xx;xx;xx
2105	ステータス 経過 ^{*1)} 、現在の CO ₂ オフセット + 待ち時間 [分:秒] 自動スタートおよび前の開始 日付と時刻	25 バイト / 十進値 フォーマット xx;xx.x;+xxx:xx;xx.xx.xx;xx:xx
2140	ステータス ガスボトル切替え 装置 CO ₂ を読み取る ^{*3)}	2 バイト / 十六進 フォーマット xx
2141	ステータス ガスボトル切替え 装置 0 ₂ を読み取る ^{*3)}	2 バイト / 十六進 フォーマット xx
2300	エラーメモリを選択 (現在のエラー) ^{*4)}	~ 241 バイト / 十六進値 フォーマット 該当の章を参照
2301	エラーメモリを選択 (過去エラー) ^{*4)}	~ 241 バイト / 十六進値 フォーマット 該当の章を参照
2400	データログに保存されている データのデータ確認 (スタート) ^{*5)}	~ 224 バイト / 十六進値 フォーマット 該当の章を参照
2401	データログに保存されている その他のデータの確認 ^{*6)}	224 バイト / 十六進値 フォーマット 該当の章を参照
2402	最近のデータログ確認の (反復) データ確認 ^{*7)}	224 バイト / 十六進値 フォーマット 該当の章を参照
2410	書込みサイクル データログを 時間:分:秒で選択	8 バイト / 十進値 フォーマッ ト xx:xx:xx

*1) 表を参照 注 ステータス プロセス 消毒および自動スタート。

*2) レベルごとに各2 バイト。

*3) ボトル A オン (0x01)、ボトル B オン (0x02)、圧 ボトル A OK (0x10)、圧ボトル B OK (0x20)。

*4) エラーメモリに関する詳細は、13.5章を参照。

*5) リードポインタを最初の入力に据える、最大7つの入力を読み取る。

*6)次の7つの入力を送る。リードポインタを自動で次の新しい入力に据える、max.7 つの入力を読み取る。

*7) もっとも最近のテレグラムの入力を、改めて送信。接続障害が起こった場合に、ご利用いただくことができます。

Bit	消毒	自動スタート
0x00	Steri-Run オフSteri-Run オフ	自動スタートオフ
0x01	初期化	初期化
0x02	開扉時間を待つ	開扉時間を待つ
0x03	閉扉を待つ	閉扉を待つ
0x04	スタート	スタート
0x05	ヒーティング	ヒーティング
0x06	ストップ	逆電流調整を行う
0x07	コンデンゼーション	待ち時間 1
0x08	冷却	公差範囲を据える
0x09	乾燥	安定した湿度を築く
0x0A	リリースを待つ	逆電流調整を行う
0x0B	中断	待ち時間 2
0x0C	-	オフセットを算出する
OxOD	-	オフセットを選択し確認する
0x0E	-	リリース
0x0F	-	中断

注*3) ステータス プロセス 消毒および自動スタート:

構造 エラーメモリ

エラーメモリには、22 つのエラー表示が保存されています。データ確認は、22 つの記録を 夫々コロンで区切った形で返答が行われます。以下のコマンドでデータ確認が可能です:

- データ確認: ?:2300:00::cc<CR> エラーメモリ内の最新入力情報 11 つを読み込む
- データ確認: ?:2301:00::cc<CR> エラーメモリ内の最初入力情報 11 つを読み込む

これらの記録は11バイトで、21 つの ASCII 記号に移行する前に暗号化されます。例えば、バイト 0x23 から ASCII 記号 0x32 (「2」)と 0x33 (「3」)のようにです)。
・バイト 1 は、1 つの記号から成っています。
・バイト 2 ~ 11 は、2 つの記号から成っています。

すなわち返答は、1+(10 x 2) = 21 バイト、プラス区切りのコロンで構成されています。 記録一つの中には、日付、時刻、エラーのあるループ、装置ステータス、エラー表示情 報が入っています。

返答の例:

!:2300:fb:10b01060f372280000002:20b01060f38100001....:80

- 記録1: !:2300:fb:10b01060f37228000002: (21 バイト)
- 記録 2: 20b01060f38100001.....:80 (記録 2 は、記録 1 の 01060 バイトおよび 区切り記号 [1 バイト] の後に続く)

パターン 記録構造 エラーメモリ:

	00	0b	01	06	0f	37	22	8000	0002
エラーを含むループのタイプ									
0x00 = 温度ループのエラー	00								
日付 / 日 0x0b = 月の11日		0b							
日付 / 月 0x01 = 1月			01						
日付 / 年 0x06 = 2006				06					
時刻 / 時 0x0f = 15時					Of				
時刻 / 分 0x37 = 55 分						37			
時刻 / 秒 0x22 = 34 秒							22		
装置ステータス 0x8000 = 装置エラー有								8000	
エラー表示 0x0002 = 実測値が高すぎる									0002

この記録内には、以下の情報も移行されます:

- ・作成日 2006 年1月11日、15:55:34 時。
- ・装置エラーがあります。実測温度 が高すぎます。

エラー表示の一覧は Hex コーデックにて

Hex ⊐ード	説明 / タイプ
0x00	ループ 温度
0x01	ループ CO ₂
0x02	ループ 0 ₂
0x07	水位
0x08	一般装置ステータス

エラー表示の一覧はBit コーデックにて 一般装置ステータス、ループ温度、CO₂:

Bit	一般装置ステータス
0x0002	装置ドアの開扉時間が長すぎ
0x0004	ディスプレイの反応がない
0x0008	パラメータ メインボードが妥当でない(EEPROM 欠陥)
0x0010	データログの欠陥(装置は引き続き作動可能)
0x0020	消毒 / Steri-Run内でのエラー
0x0040	Steri-Run 中の電気供給がない
0x0080	自動スタート内でのエラー
0x0100	ADC テスト失敗
0x0400	ファンエラー
0x1000	IR センサー交換(Info)
0x2000	自動スタートオン(Info)
0x4000	消毒 オン(Info)
0x8000	装置エラーあり(Info)
Bit	エラーステータス ループ温度
0x0001	センサー破損
0x0002	実測値 大きい
0x0004	実測値 小さい
0x0008	実測値 妥当ではない
0x0010	校正値 大き/小さすぎ
Rit.	エラーステータス ループ 00。

Bit	エラーステータス ループ CO ₂
0x0001	センサー破損
0x0002	実測値 大きい

Bit	エラーステータス ループ CO ₂
0x0004	実測値 小さい
0x0010	校正値 大き/小さすぎ
0x0020	接続障害(センサー)
0x0040	接続障害(ガスボトル切替え装置)
0x0080	ガスなし、ボトル A と B が空
0x0200	ガスボトル A 空
0x0400	ガスボトル B 空

ループ 02 および水位:

Bit	エラーステータス ループ O2
0x0001	センサー破損
0x0002	実測値 大きい
0x0004	実測値 小さい
0x0020	接続障害(センサー)
0x0040	ガスボトル切替え装置 の反応がない
0x0080	ガスなし、ボトル A と B が空
0x0200	ガスボトル A 空
0x0400	ガスボトル B 空
Bit	エラーステータス 水位
0x0001	水なし

データログ構造

データログが 10.000 個までの入力情報を保存します。記録インターバルの設定によって は(秒単位)、例えば 10,000 秒設定(デフォルト)の場合、約5日間の記録を残すこと が可能です。

データログ内には、以下の情報が保存されます:

· 重要なユーザーアクション、システムイベント、エラー表示情報。

・インキュベーション運転中の3ループでの計測データ

データログは、以下のコマンドにてデータ確認ができます:
 データ確認: ?:2400:00::cc<CR>
 リードポインタをデータログのリードポインタを最も古い入力情報と最初の記録アウトプットに据える。

 データ確認: ?:2401:00::cc<CR>
 以下の記録のアウトプット、リードポインタは自動で少しずつ過去から最新の入力情報へと移動します。

データ確認: ?:2402:00::cc<CR> 最近読まれたデータの再アウトプット、リードポインタはこの コマンド時には動きません。このコマンドで、 接続エラー後のデータ損失を防ぐことができます。

データ確認コマンドはそれぞれ、区切り記号なしで、7 つまでの記録を使い返答されま す。これらの記録は 16 バイトで、32 つの ASCII 記号に移行する前に暗号化されます。 すなわち、例えばバイト 0x23 から、ASCII 記号: $0x32([2_]) \ge 0x33([3_])$ 。 よって、返答 1 つは 7 × 16 = 112 バイトまで、すなわち 224 の ASCII 記号から構成さ れます。 記録一つには常に、日付、(秒表記なし)、装置ステータス、データログ入力情報のタイ プが含まれます(バイト 0 ~ 7 または ASCII 記号 0 ~ 15)。 入力情報によって、ループの実測または設定値、もしくはその他のパラメータを入寮す ることが可能です(バイト 8 ~ 15、または ASCII 記号 16 ~ 31)。

返答の例:

!:2400:e0:010b01060f3700000177002800d40000110b01060f3800000172003200d20352... ...:80

- 記録1 !:2400:e0:010b01060f3700000177002800d4000011
 - (32 バイト ASCII 記号)
- 記録 20b01060f3800000172003200d20352...80(記録 2 は、記録 1 の 32 バイトの後に続く)

パターン 記録構造 データログ:

データログの記録タイプ	01	0b	01	06	Of	37	0000	0177	0028	00d4	0000
0x01 = 60 秒毎に保存される標準記録	01										
日付 / 日 0x0b = 月の11日		0b									
日付 / 月 <u>0x01 = 1月</u>			01								
日付 / 年 0x06 = 2006 				06							
時刻 / 時 0x0f = 15時					0f						
時刻 / 分 0x37 = 55 分						37					
装置ステータス 0x0000 = 問題なし							0000				
タイプ別、 現在の温度 摂氏1/10℃単位 0x0177 = 37.5 ℃								0177			
タイプ別、 現在のCO2含有量 1/10%単位 0x0028 = 4.0 % のCO ₂									0028		
タイプ別、											
現在のO 含有量 1/10%単位 0x00d4 = 21 2 %の O ₂											
<u>rH用</u>										00d4	
1/10%単位 常に 0x0000 = 0%											0000

この記録内には、以下の情報も移行されます:

- · 作成日 2006 年 1 月 11 日、15:55 時。
- ・装置ステータスは特に異常を表示していません。
- ・温度 は 37.5 °C。
- ・ガス濃度 4.0% CO₂、21.2% O₂。

留意コード例: コード例は、本章の接続の説明部分にございます。

イベント入力情報の一覧は Bit コーデックにて イベント入力情報一覧パート I:

コード	イベント	特記事項(バイト 8 ~ 15)
0x01	全ループの設定値 (周期的、分サイクル)	現在值 温度、CO ₂ 、O ₂ 、rH
0x02	設定値変更 (新しい段階を始める際に)	設定值 温度、CO ₂ 、O ₂ 、rH
0x10	設定値 温度 変更する	設定值 温度、CO ₂ 、O ₂ 、rH
0x11	設定値 CO ₂ 変更する	設定值 温度、CO ₂ 、O ₂ 、rH
0x12	設定値 0 ₂ 変更する	設定值 温度、CO ₂ 、O ₂ 、rH
0x20	新エラー 温度	ステータス / エラー登録 温 度、CO ₂ 、O ₂ 、rH
0x21	新エラー CO ₂	ステータス / エラー登録 温 度、CO ₂ 、O ₂ 、rH
0x22	新エラー 0 ₂	ステータス / エラー登録 温 度、CO ₂ 、O ₂ 、rH
0x2F	新エラー System	ステータス / エラー登録 温 度、CO ₂ 、O ₂ 、rH
0x30	ネットワークリセット	設定值 温度、CO ₂ 、O ₂ 、rH
0x31	カバーが開いています	現在の実測値 温度、CO ₂ 、O ₂ 、 rH
0x32	扉が閉じている	現在の実測値 温度、CO ₂ 、O ₂ 、 rH
0x40	顧客校正 温度	校正レベル(2 バイト)、温度 旧、温度 新(各2 バイト)
0x41	顧客校正 CO2	校正レベル(2 バイト)、CO ₂ 値 旧、CO ₂ 値 新(各2 バイト)
0x42	顧客校正 02	校正レベル(2 バイト)、0 ₂ 値 旧、0 ₂ 値 新(各2 バイト)
0x50	自動スタート Start	ステータス / エラー登録 温 度、CO ₂ 、O ₂ 、rH
0x51	自動スタート完了	現在の実測値 温度、CO ₂ 、O ₂ 、 rH

イベント入力情報一覧パート II:

コード	イベント	特記事項(バイト 8 ~ 15)
0x52	自動スタート エラー発生にて 終了	ステータス / エラー登録 温度、CO ₂ 、O ₂ 、 rH
0x53	自動スタート手動停止	ステータス / エラー登録 温度、CO ₂ 、O ₂ 、 rH
0x60	Start Steri-Run	ステータス / エラー登録 温度、CO ₂ 、O ₂ 、 rH
0x61	Steri-Run 完了	現在の実測値 温度、CO ₂ 、O ₂ 、rH
0x62	Steri-Run エラー発生にて終了	ステータス / エラー登録 温度、CO ₂ 、O ₂ 、 rH
0x63	Steri-Run 手動停止	ステータス / エラー登録 温度、CO ₂ 、O ₂ 、 rH
0x70	ガスモニター ボトル A 空	ステータス ガスモニタリング (2 バイト)、 4 バイト 空
0x71	ガスモニター ボトル B 空	ステータス ガスモニタリング (2 バイト)、 4 バイト 空
0x72	ガスモニター 手動切替え	ステータス ガスモニタリング (2 バイト)、 4 バイト 空
0x90	低湿度 Start	現在の実測値 温度、CO ₂ 、O ₂ 、rH
0x91	低湿度 Stop	現在の実測値 温度、CO ₂ 、O ₂ 、rH
0xe0	データログ消去	現在の実測値 温度、CO ₂ 、O ₂ 、rH
0xff	データログへの最新の入力情報	情報なし。日付、時間、ステータスに関す る情報もなし

コード例 データログ

データログの入力情報は、1つ16 バイトの大きさで、以下の通りの構成となっています:

- 1. バイト: イベント表示 (例えば、扉が開いている 0x31、測定値入力 0x01)
- 2. バイト:入力日
- 3. バイト:月
- 4. バイト:年
- 5. バイト:時間
- 6. バイト:分
- 7. ~ 8. バイト:装置ステータス
- 9. ~ 16. バイト: イベントに関する様々なデータ

データログデータ確認の機能

以下のデータログ選択におけるコード例では、6つの機能が使用されます:

- ahex
 - // 受信した ASCII 記号を十六進数に変換、
- send_telegramm
 // データログに問い合わせを送信、
- get_telegramm
 // データログより返答を受信、
- time_2_str
 // 十六進値から時刻フォーマットの ASCII 記号を作成、
 num_2_string
 - // 十六進値からファイルに入力するために ASCII 記号を作成、
- read_datalogger
 // 受信データを処理しファイルに記入。

データログデータ確認のコード例

char ahex (char a)

```
char ahex(char a)
{
    char i;
    char hexa[16]= 「0123456789abcdef」;
    for (i = 0; i < 16; i++)
        if (a == hexa[i])
        return (i);
    return 0;
}</pre>
```

send_telegramm

```
void send_telegramm(char *p)
   ł
  char string [15];
  unsigned char bcc = 0xFF;
  char i;
// テレグラムを一緒にコピー
  strncpy (&string[0], [?:xxxx:00::00\r], 14);
// 4桁のアドレスを挿入
  strncpy (&string[2], p, 4);
// チェック数を算出: 全バイトの逆転 XOR
// チェック数なし、そして <CR>
  for (i = 0; i < 11; i ++)
     bcc = (bcc^string[i]);
// チェック数 をコピー
  string[11] = hexa(bcc/16);
  string[12] = hexa(bcc%16);
```

// テレグラム送信 ComWrt (COM_NR, string, 14); return; }

get_telegramm

```
int get_telegramm(char *p)
{
    int reading_count = 0;
// テレグラムの記号による選択
    do
        ComRd(COM_NR, &p[reading_count], 1);
// 受け取るまで <CR>
    while ((p[reading_count++] != '\r'));
// 返却 = 受信した記号数
    return (reading_count);
```

time_2_str

```
char time_2_str (int z, char * b)
{
     char i;
// 数字を二つ出力
     for (i = 1; i >= 0; i--) {
     // 値を算出
        b[i] = z%10+0x30;
     // 基準値を下げる
     z = z/10;
     }
     return (2);
     }
```

num_2_string

```
rest = z % 10;
// 小数点以下を切り落とす
  z = z / 10;
// 小数点以上の数字を算出しコピー
  for (i = 0; i < 12; i++)
// 値を算出
  a[i] = z%10+0x30;
// 基準値を下げる
  z = z/10;
// 数字を完全にコピーしましたか?
  if (z == 0) break;
  }
  for (; i \ge 0; i-)b[l++] = a[i];
// コンマ以下の数を算出しコピー
  b[l++] = ',';
// 値を算出
  b[l++] = rest%10+0x30;
  return (I);
  ł
```

read_datalogger

```
int read_datalogger ()
ł
#define SIZE_DATA2 16
#define EVENT_STATUS 0x01
unsigned char buffer[300], string [300];
unsigned char zahlenstring [150], datestring, timestring;
unsigned char excelstring [150];
unsigned char len, h, i;
unsigned int read_count, status;
#define EVENT_DATA. END 0xFF
char data;
int GetTele = 0
GetError = 0,
// ファイルのタイトル行に記載
WriteFile (FileHandle, 「Date; Time; Comment; Temp Act.; CO2 Act.; O2
Act.;rH Act.;Temp Set;CO2 Set;O2 Set;rH Set;\nj , 85);
// エンドロスループ
while (1)
// データログを初めに据えて読みとる
   if (!GetTele) {
   send_telegramm (「2400」);
   }
   else{
// 更に記録を読み取る
   send_telegramm (「2401」);
   ł
   len = get_telegramm (buffer);
// テレグラムが受信できません
```

```
if (!len) {
      GetError ++;
// 新しい問合せ
   send_telegramm (「2402」);
   len = get_telegramm (buffer);
// またしてもテレグラムが受信できません
   if (!len) return 1;
   }
// テレグラムカウンターを上げる
   GetTele ++;
// 送付した使用データの長さ
   len = (ahex(buffer[7]) * 0x10 + ahex(buffer[8])) / 2;
// ASCII ストリングを利用可能な数字のストリングに変換
   for (i = 0; i < (string); i++)
   zahlenstring [i] = (ahex(buffer[10 + (2*i)]) * 0x10 +
   ahex(buffer[11 + (2*i)]));
// 送付されたデータパッケージの算出
   data = ((len) / SIZE_DATA2);
// 全データパッケージの評価
   for (i = 0; i < data; i++)9{
   len = 0:
// 時刻と日付をファイルに記載
   len += time_2_str (zahlenstring[1+i*SIZE_DATA2],
   &excelstring[len]);
   excelstring[len ++] = '.' ;
   len += time_2_str (zahlenstring[2+i*SIZE_DATA2],
   &excelstring[len]);
   excelstring[len ++] = '.' ;
   len += time_2_str (zahlenstring[3+i*SIZE_DATA2],
   &excelstring[len]);
   excelstring[len ++] = ';' ;
   len += time_2_str (zahlenstring[4+i*SIZE_DATA2],
   &excelstring[len]);
   excelstring[len ++] = ':' ;
   len += time_2_str (zahlenstring[5+i*SIZE_DATA2],
   &excelstring[len]);
   excelstring[len ++] = ':' ;
   len += time_2_str (0, &excelstring[len]);
   excelstring[len ++] = ';';
   switch (zahlenstring[i*SIZE_DATA2]) {
      case EVENT_STATUS:
// サイクルの入力情報を確認 装置エラー
   status = zahlenstring[6+i*SIZE_DATA2]*0x100+
   zahlenstring[7+i*SIZE_DATA2];
      if (status & INFO_ERROR) {
         str_cpy (&excelstring[len], 「Error active;」, 13);
         len += 13;
      }
      else{
```

```
// 全装置エラー データ確認(参照 「イベント入力情報の一覧はBitコーデックにて
イベント入力情報一覧パート I:」上 頁 12-14)
         if (status & DOOR_LONG) {
            str_cpy (&excelstring[len], 「Door open too long;」,
            19);
            len += 19;
         else {
            if (status & DOOR_OPEN) {
               str_cpy (&excelstring[len], 「Door open;」, 10);
               len += 10;
               }
            }
// 残りの装置エラー データ確認
//
//
//
//
//
// そして最後にサイクルの実測値入力 装置エラーなしで
// データ確認
else{
         str_cpy (&string[string], 「ok;」, 3);
         string += 3;
      }
   }
// 数字ストリングの実測値をエクセルストリングにコピー
   len += num_2_str ((zahlenstring[8+i*SIZE_DATA2]*0x100+
   zahlenstring[9+i*SIZE_DATA2]), &excelstring[len]);
   excelstring[len ++] = ';' ;
   len += num_2_str ((zahlenstring[10+i*SIZE_DATA2]*0x100+
   zahlenstring[11+i*SIZE_DATA2]), &excelstring[len]);
   excelstring[len ++] = ';' ;
   len += num 2 str ((zahlenstring[12+i*SIZE DATA2]*0x100+
   zahlenstring[13+i*SIZE_DATA2]), &excelstring[len]);
   excelstring[len ++] = ';' ;
   len += num_2_str ((zahlenstring[14+i*SIZE_DATA2]*0x100+
   zahlenstring[15+i*SIZE DATA2]), &excelstring[len]);
   excelstring[len ++] = ';' ;
// これ以降設定値を入力
   len += num_2_str (SollTemp, &excelstring[len]);
   excelstring[len ++] = ';' ;
   len += num 2 str (SollCO2, &excelstring[len]);
   excelstring[len ++] = ';'
   len += num_2_str (SollO2, &excelstring[len]);
   excelstring[len ++] = ';' ;
   len += num_2_str (SollrH, &excelstring[len]);
   excelstring[len ++] = ';' ;
   excelstring[len] = '¥n' ;
   len += 1;
   WriteFile (FileHandle, excelstring, len);
```

```
break;
// これ以降残りのイベントのデータ確認
   case EVENT_FORMAT_DATALOG:
      WriteFile (FileHandle, excelstring, len);
      WriteFile (FileHandle, 「Data logger erased;\n」, 20);
      break;
   case EVENT POWER ON:
// 設定値のアップデート
      SetTemp = zahlenstring[8+i*SIZE_DATA2]*0x100+
      zahlenstring[9+i*SIZE_DATA2];
      SollCO2 = zahlenstring[10+i*SIZE DATA2]*0x100+
      zahlenstring[11+i*SIZE_DATA2];
      Soll02 = zahlenstring[12+i*SIZE_DATA2]*0x100+
      zahlenstring[13+i*SIZE_DATA2];
      SollrH = zahlenstring[14+i*SIZE_DATA2]*0x100+
      zahlenstring[15+i*SIZE_DATA2];
      WriteFile (FileHandle, excelstring, len);
      WriteFile (FileHandle, 「Power on;\n」, 10);
      break;
   case. .
// ここで全イベントのデータ確認(参照 「イベント入力情報の一覧は Bit コーデック
にて イベント入力情報一覧パート I:」上 頁 12-14)
// 中断 0xFF とあればデータログの最後です
   case 0xFF:
  WriteFile (FileHandle, [End;\n], 5);
   ł
}
return 0;
ł
```

プログラム STERI-CYCLE i160 LK/ STERI-CYCLE i250 LK

このプログラムは、装置と接続されたパソコン間におけるデータコミュニケーションの 処理のため、ユーザーインターフェース (英語のメニュー名のみ)を設けています。



このプログラムは、下記の用途で使用されます:

- エラー表示情報(エラーログ)の選択とアーカイブ。記録情報は、メタフォーマット *. CSV で保存されます。
- ・イベント入力情報(データログ)の選択とアーカイブ。記録情報は、メタフォーマット *. CSV で保存されます。
- Thermo Fisher Scientific テクニカルサービスへ送付される、サービスファイルの 作成。サービスファイルの情報を基に、体系的なエラー調査を行う事が可能です。
 記録は独自フォーマット *. SRF として保存されます:

STERI-CYCLE i160 LK & STERI-CYCLE i250 LK のインストール

- 1。インストールをスタート:
- · CD-ROM 上のサブディレクトリ PROGRAMS die Datei SETUP.EXE をダブルクリックで開らきます。

	bin
	license
	supportfiles
	dp.pmf
	nidist.id
-	setup.exe
	setup.ini

2。インストールリストをプログラム用に定めます。

Destination Directory Select the primary installation directory.	
All software will be installed in the following locations. To install so different location, click the Browse button and select another direr	ftware into a ctory.
Target directory for application C:\Program Files (x86)\TFS\	Browse
T and dealer for Mating II also and a financial	
Target directory for National Instruments software	Browse
C:\Program Files (x86)\National Instruments Sottware	

- 3。記載されているインストール手順の通りに:
 - ・ライセンス規定に同意する、
 - ・インストールの範囲を決める、
 - インストール完了の表示が出たら、インストールインターフェースを閉じ、コン ピューターを再起動します。

STERI-CYCLE i160 LK & STERI-CYCLE i250 LK の使用

ユーザーメニューの構造:

- ユーザーインターフェースは、主要メニュー2つから構成されています:
- ・ 両機能が付いた MAIN:
 - ・プログラムバージョンの表示: FIRMWARE VERSION
 - ・プログラム終了のボタン:QUIT

Thermo Fisher CO2 Incubators COM: NoCOM	_ _ X
MAIN	GENERAL
50135297 V2.7 20.05.2014 Device Name no com.	Thermo Fisher SCIENTIFIC
Date: no com. Firmware version: no com	Quit

・サブメニューの付いた GENERAL:

- · 伝達速度の設定やシリーズ接続の選択をする PRESETTING、
- · パソコンからインキュベータへのコミュニケーション接続をテストする TEST COM、
- · ご希望のタイムゾーンにおける日付と時間の設定を行う DATE & TIME、
- ・エラー表示の選択をする ERROR LOGGER、
- · イベント入力情報の選択をする DATA LOGGER、
- ・エラー情報の選択とサービスファイルの作成を行う SERVICEFILE
- · インキュベータの装置パラメータへのアクセスを管理する PASSWORD

機能とユーザーメニュー:

PRESETTING

サブメニュー PRESETTING では、伝達速度の設定やシリーズ接続の選択を行うことができます。

Thermo Fisher CO2 Incubate	ors COM: NoCOM				
	MAIN			GENERAL	
PRESETTING	TEST COM	DATE & TIME	LOGGER	SERVICEFILE	PASSWORD
RESELTING Serial Baudrate 57 600 (chang save setting	rESI COM presetting: Com-Number com-Number e ato file	DATE & TIME	LOGGER	SERVICEFILE	PASSWORD

- 1。9600~115200 ボーの間で伝達速度を選択。
- パソコンのシリーズ接続を選択。USB ドライバがインストールされている場合は、 USB 接続端子に分配されている(仮) Com ポートを選択することができます(参照 「USB インターフェース」上 頁 12-1)。
- 3。設定を引き継ぐ:
 CHANGE キーを押します。
- 4。設定を保存します (Ini ファイル): · SAVE TO FILE キーを押します。

留意 伝達速度:

伝達速度の設定は、ユーザーメニュー PRESETTING と装置において同様の値である必要 があります!

TEST COM

サブメニュー TEST COM では、サブメニュー PRESETTING にて設定した値のコミュニケー ション接続をテストすることができます。

	MAIN			GENERAL	
PRESETTING	TEST COM	DATE & TIME	LOGGER	SERVICEFILE	PASSWORE
command (!) / req. 1 2 2 address (e.g. 20 2010 attachment (e.g. string sent string sent string received	nest (?) 110) . +037.00;)				send
]

- 1。インキュベータの測定可能な最新温度値のデータ確認例:
- ・データ確認:?(プリ設定、変更不可)
- ・ アドレス: 2010(アドレス 温度値:
- 2。インキュベータヘデータ確認を送信:
- · SEND キーを押します。
- インキュベータが返答ストリングを送り返して来たら、インキュベータへの接続が されているという事になります。
- ・ 接続ができない場合は、エラーダイアログが表示されます:



- ・3. エラーダイアログを閉じる:
- · OK キーを押します。

DATE & TIME サブメニュー DATE & TIME では、日付と時間をご希望のタイムゾーンに合わせることがで きます。

mo Fisher CO2 Incubators COM: NoCOM				
MAIN			GENERAL	
PRESETTING TEST COM	DATE & TIME	LOGGER	SERVICEFILE	PASSWORD
device date:				
(dd.mm.yy) device time:				
(hh.mm:ss)				
(dd.mm.yy hh.mm.ss)				

- 1。両方とも、TT.MM.JJ(日、月、年)の順で入力してください。
- 2。入力値を適用させる:
- · SET キーを押します。

ERROR LOGGER

サブメニュー ERROR LOGGER では、ユーザーインターフェースのテキストボックスにエ ラー表示を読み込ませることができます。

Thermo Fisher CO2 Incubators CO	M: NoCOM				_ _ ×
	MAIN			GENERAL	
PRESETTING	TEST COM	DATE & TIME	LOGGER	SERVICEFILE	PASSWORD
DATA LOGGER	ERROR LOGGER				
	date	time	Control loop	Loop n	nessage
read Amount data 0 Amount telegrams					
Amount errors					save to file

記録情報は、メタフォーマット *. CSV にて保存することができます。 ・記録情報をファイルとして保存する:

· SAVE TO FILE キーを押します。

DATA LOGGER

サブメニュー DATA LOGGER では、ユーザーインターフェースのテキストボックスにエ ラー表示を読み込ませることができます。

記録情報は、メタフォーマット *. CSV にて保存することができます。

- 1。記録を読み込む:
- · READ キーを押します。
- 2。データ移行の経過は、3つのテキストボックスにて表示がされます:
- · AMOUNT DATA: データ移行される記録の総数
- · AMOUNT TELEGRAMS: 移行されるテレグラムの数
- · AMOUNT ERRORS: 送信されたエラーメッセージの数。

留意 データ移行にかかる時間:

データログは最大1万個の記録情報を有することができるということがあり、パソコンへのデータ移行には時間がかかる可能性があります。

SERVICEFILE

サブメニュー SERVICEFILE で、インキュベーターのエラー情報の読み込みとサービス データの作成ができます。サービスデータは、独自フォーマットの *. srf にて保存され ます。サービスデータは、エラー分析の為 Thermo Fisher Scientific 社のテクニカル サービスに転送されます。

	MAIN					GENERAL	
PRESETTING	TEST COM	DATE & 1	TIME	LOGGE	ER S	SERVICEFILE	PASSWORE
	Öffnen					×	
	Directory History: C:\					•	
	Suchen in:	SDisk (C:)		•	+ 🗈 💣 🗊 🕶		
	œ.	Name	^		Änderungsdatum	Тур	
	~	Benutzer			21.02.2014 13:24	Dateio	
	Zuletzt besucht	CDs			02.04.2014 11:19	Dateio	
create servicefile		📕 dell			04.04.2011 20:36	Dateio	
	Desktop	b Doku			11.06.2014 13:54	Dateio	
Amount data	(Second p	ownload 🔒			24.01.2014 11:54	Dateio	
0	-	DPSEDriver			14.11.2013 14:25	Dateio	
	Bibliotheken	drivers			08.01.2014 07:13	Dateio	
Amount telegrams		FILES			22.03.2012 13:17	Dateio	
0		Hotfix			10.02.2012 00:14	Dateio	
	Computer	🗼 Intel			17.01.2011 08:44	Dateio	
		📕 Kpcms			16.03.2012 08:01	Dateio	
	Network	DM-DELSB			28.09.2012 07:20	Dateio	
	INELZWEIK	Perfl ons			14 07 2009 05-20	Datein *	
		•	m			•	
		Dateiname: serv	ice.srf		•	ОК	
		Dateityp: (*.si	rf)		•	Abbrechen	
		_	_	-	_		

- 1。サービスファイルを作成する:
 - · CREATE SERVICEFILE キーを押します。
- · Windows のダイアログボックスに、保存のためファイル名とリストを定めます。
- 2。保存プロセスを開始:
 OK キーを押します。

留意作成にかかる時間: 装置情報の収集とサービスデータの作成には、時間がかかる可能性があります。

<mark>12 添付物 データ通信</mark> プログラム STERI-CYCLE i160 LK / STERI-CYCLE i250 LK

PASSWORD

サブメニュー PASSWORD は、Thermo Fisher Scientificのサービス職員のみ使うことができます。

mo Fisher CO2 Incubate	ors COM: NoCOM				
	MAIN			GENERAL	
PRESETTING	TEST COM	DATE & TIME	LOGGER	SERVICEFILE	PASSWORD
		PASSWORD			
			ENTER		

13-

装置ブック

既にした作業を、ここに入力してください:

装置タイプ:		注文番号:		
工場番号:		サービス番号:		
設置場所:		オペレーターの注:		
行った作業:	注:	·	日付	署名

13 装置ブック

Thermo Scientific のコンタクトデータ

Thermo Fisher 社の世界中の営業組織

住所 ドイツ: Thermo Electron LED GmbH Robert-Bosch-Straße 1 D - 63505 Langenselbold

ドイツからのお問い合わせ:

Tel 営業 0800 1 536376 Tel サービス 0800 1 112110 Fax 営業・サービス 0800 1 112114 E-Mail info.labequipment.de@thermofisher.com ヘルプデスクEメール: service.lpg.germany.de@ThermoFisher.com

Enquiries from Europe, Middle East and Africa:

Tel. + 49 (0) 6184 / 90-6940 Fax: + 49 (0) 6184 / 90-7474 E-Mail info.labequipment.de@thermofisher.com

Postal address USA: Thermo Scientific 275 Aiken Road Asheville, NC 28804 USA

Enquiries from North America: Phone +1 800-879 7767 +1 800-879 7767 Fax +1 828-658 0363 Email: info.labequipment@thermofisher.com

Enquiries from Latin America: Phone +1 828-658 2711 Fax +1 828-645 9466 Email: info.labequipment@thermofisher.com

Enquiries from Asia Pacific: Phone +852-2711 3910 Fax +852-2711 3858 Email: info.labequipment@thermofisher.com

Enquiries al address USA:

Thermo Scientific 275 Aiken Road Asheville, NC 28804 USA

Enquiries from USA/Canada

Sales: +1 866 984 3766 Service+1 800 438 4851

Enquiries from Latin America Sales: +1 866 984 3766 Service:+1 866 984 3766

Enquiries from Asia: China Sales: +86 10 8419 3588 Service:Toll free 8008105118 Support Mobile 4006505118 or +86 10 8419 3588

India

Sales: +91 22 6716 2200

Service: Toll free 1 800 22 8374 or +91 22 6716 2200

Japan Sales: +81 45 453 9220 Service:+81 45 453 9224

Enquiries from the Rest of Asia/Australia/New Zealand Sales: +852 2885 4613 Service:+65 6872 9720

Enquiries from Countries not listed / Rest of EMEA Sales: +49 6184 90 6940 or +33 2 2803 2000 Service:+49 6184 90 6940

Enquiries from Europe: Austria Sales: +43 1 801 40 0 Service:+43 1 801 40 0

Belgium Sales: +32 53 73 4241 Service:+32 53 73 4241

Finland/Nordic/Baltic countries Sales: +358 9 329 100 Service:+358 9 329 100

France

Sales: +33 2 2803 2180 **Service:**+33 825 800 119

Germany: Postal address Germany: Thermo Electron LED GmbH Robert-Bosch-Straße 1 D - 63505 Langenselbold Phone Sales Toll free 0800 1 536 376 or +49 6184 90 6940 Service Toll free 0800 1 112110 or +49 6184 90 6940 E-Mail info. labequipment. de@thermofisher. com Italv Sales +39 02 95059 341 Service+39 02 95059 250 Nether lands Sales +31 76 579 5555 Service+31 76 579 5639 Russia/CIS Sales +7 812 703 4215 Service+7 812 703 4215 Spain/Portugal Sales +34 93 223 0918 Service+34 93 223 0918 Switzerland Sales +41 61 716 7755 Service+41 61 716 7755 UK/Ireland Service+44 870 609 9203 Sales +44 870 609 9203

[®] 2013 Thermo Fisher Scientific Inc. 全権利を留保します 本書記載のすべての商標は Thermo Fisher Scientific 社およびその関連会社の独占的財産です。表示の仕様、条 件、価格に、拘束力はありません。国によっては、お取り扱いのない製品もございま す。より詳細な情報はご請求に応じてお客様現地の提携会社にて入手可能です。 14 Thermo Scientificのコンタクトデータ

A

Cell locker 用 6 ドア気密スクリーン

Cell locker 用 6 ドア気密スクリーンは Cell locker 装置用に設計された 6 室に仕切られ た内扉です。

Cell locker 用6ドア気密スクリーンを使用すると、CO₂インキュベーターの内部を6つ のサンプル室(Cell locker)に分割することができます。サンプルの取り出しができる ように、前面プレートにはガラスドア(通路開口部)が設けられており、各ドアが開閉 できるようになっています。



Cell locker 用 6 ドア気密スクリーン



CELL LOCKER

留意 インキュベーターの性能の変化 6ドア気密スクリーンの設計上、インキュベーターの性能が変化します(仕様を参照)。 A Cell locker用6ドア気密スクリーン 棚板の設置

棚板の設置



棚受け(1)をサポートレールの下部、中央部、上部にある長方形の穴に挿入します。 上部および中央部の位置に長い(一本につながった)ガイドレール付きの有孔棚板を設 置します。



図: 長い(一本につながった)ガイドレール付きの有孔棚板

シリコンストッパー(1)を取り外さないでください。

ガイドレール(2)を取り付けてから、長いガイドレール付きの有孔棚板を設置してくだ さい。

長いガイドレールは必要に応じて取り外すことができます。

中間がない2分割されたガイドレール(1)付きの有孔棚板を下部の位置に設置します。



図: 下部に配置する2分割ガイドレール付きの有孔棚板

2分割ガイドレールは取り外すことができません。



図: 有孔棚板の設置

有孔棚板をレールの後方にスナップして取り付けます。有孔棚板をレールの後方にス ナップして取り付けます。ドアを閉じるためには、棚板の底面にある金属製のタブを レールの後ろに固定する必要があります。

6ドア気密スクリーンを閉じます。

A Cell locker用6ドア気密スクリーン 棚板の設置

取扱説明書

Cell Locker 50156300 リビジョン D 2021 年 7 月



前書き

© 2021 Thermo Fisher Scientific Inc. 全権利を留保します。

Thermo Fisher Scientific社は、製品の操作に使用する目的でご購入いただいたお客さまに 本書を提供しております。本書は著作権によって保護されており、Thermo Fisher Scientific 社の書面による許可がない限り、本書の全部または一部を複製することは固く禁じられていま す。

本書の内容は予告なしに変更されることがあります。

Thermo Fisher Scientific社は、本書が完結していることや正確で誤りのないことを明言す るものでもなければ、一切の責任を負うものでもありません。また、本書の内容に忠実に従っ たとしても、本書の使用に起因するいかなる誤り、脱落、損害、損失に対して法的責任を負わ ないものとします。

本書は、Thermo Fisher Scientific 社とご購入者さまとの販売契約の一部ではありません。 本書はいかなる契約条件をも規定または変更するものではありません。

本取扱説明書のドイツ語版と翻訳版とに不一致がある場合は、ドイツ語版が効力を有するもの とします。

登録商標

CELL LOCKER® は Thermo Scientific 社の登録商標です。 Thermo Scientific は Thermo Fisher Scientific 社のブランドです。

その他すべての本取扱説明書に記載されている商標は、各製造業者の独占的財産です。

Thermo Electron LED GmbH Robert-Bosch-Straße 1 D - 63505 Langenselbold ドイツ

Thermo Electron LED GmbH は以下の企業の関連会社です。

Thermo Fisher Scientific Inc. 168 Third Avenue Waltham, MA 02451 USA
コンテンツ

第	1章 1.1. 1.2. 1.3. 1.4. 1.5. 1.5. 1.5.	一般的な注意 2 安全に関する情報と記号の説明 2 安全上の注意 3 安全に使用するためのルール 3 製品保証 3 使用目的 4 1. 正しい使用法 4 2. 誤った使用法 4
第	2 章 2.1. 2.2.	機器の納入 納入品目
第	3 章 3. 1. 3. 2. 3. 3. 3. 4.	機器の詳細
第	4 章 4.1. 4.2. 4.3.	クリーニングと消毒 10 クリーニング 10 10 払拭消毒または噴霧消毒 11 11 オートクレーブ 11 12
第	5 章 5.1. 5.2.	アフターサービス 13 メンブレンフィルターの交換13 カバーガスケットの交換15
第	6 章	予備部品および摩耗部品 16
第	7 章	廃棄 17
第	8章	仕様

1. 一般的な注意

1.1 安全に関する情報と記号の説明



警告 注意しないと、死亡または重傷を負う恐れがある危険な状況を示します。



注意 注意しないと、中軽傷を負う恐れがある危険な状況を示します。

注 注意しないと、機器や周辺にある物などに損害を与える恐れがある状況を示します。

本取扱説明書で安全上の注意がある章には、このアイコンが章のタイトルの下に表示されてい ます。アイコンが機器に表示されている場合は、取扱説明書または付随する文書に記載されて いる情報に特に注意する必要があることを示しています。
アプリケーションのヒントや役立つ情報を示しています
有害液体!
火災の危険!
有害物質!
保護手袋の着用!
セーフティゴーグルの着用!
呼吸用保護具の着用!

1.2 安全上の注意

CELL LOCKER の使用は指導を受けた職員に限ります。

本取扱説明書、適応する安全データシート、衛生ガイドライン、およびオペレーターが発行し た対応する技術上の規則を使用して、対象とする機器を使用する職員向けに文書化された手順 を作成してください。作成する手順は以下の通りです。

- · CELL LOCKER 本体と使用する付属品に対する除染対策
- 特定の薬剤を処理する際に守るべき安全対策
- ・微生物学的および生物学的サンプルなどの取り扱う際の保護具の着用
- 事故の際に行う措置
- ・ 廃棄の際の乾燥時間(例えば輸送時には2時間)
- · GLP(優良試験所規範、Good Laboratory Practice)に従って作業します

1.3 安全に使用するためのルール

サンプルの重量制限に従ってください。特に総重量と棚重量の制限に注意してください。

ドアが適切に密封されていること、破損してないことなどを6か月ごとに確認してください。

有害化学物質を含むサンプルは使用しないでください。有害化学物質が欠陥のあるシールを通 して大気中に放出されたり、腐食などの欠陥を引き起こしたりする恐れがあります。

1.4 製品保証

Thermo Fisher Scientificは、以下の条件を満たした場合に限り、*CELL LOCKER* の操作上の安全 性および機能に対して1年間保証いたします。

- · CELL LOCKER が使用目的に従って、これらの操作説明書の記載どおりに操作され、技術サービスを受けている場合
- ・ 機器に修正または改良を加えていない場合
- Thermo Fisher Scientific 承認済の原初スペアパーツと付属品のみが使用されている場合 (Thermo Fisher Scientific 非承認の第三者が製造したスペアが使用されている場合は、この限定的保証は無効になります)
- 検査およびメンテナンスが指定された間隔で実行されている場合
- ・修理作業後に毎回、動作確認試験が実行されている場合

保証は CELL LOCKER の納入日から有効になります。

1.5 使用目的

1.5.1. 正しい使用法

CELL LOCKER は実験・検査に使用する装置で、特に研究用途で細胞の準備や培養、組織培養に使用されます。

高速アスピレーターを使用した場合、CELL LOCKER は 121°C (250°F) で 15 分までのオート クレーブが可能です。

1.5.2. 誤った使用法

以下のような組織、物質、液体を使用しないでください。

- ・引火性または爆発性が高いもの
- ・ 可燃性または爆発性の蒸気や粉塵を放出するもの
- ・ 毒素を放出するもの
- ・粉塵を放出するもの
- ・発熱反応を起こすもの
- 花火に含まれる物質

内面の底やユニット内部の集水槽に液体を注ぎ込まないでください。

CELL LOCKER は医療機器として使用しないでください。

2. 機器の納入

2.1 納入品目

。 説明	個数
Cell Locker	1
メンブレンフィルター	2
カバー付き作業トレイ	1
消耗品キット	1
取扱説明書	1
安全の手引き要約	1

2.2 検収

CELL LOCKER の納品時に、直ちに次の事柄を確認してください。

- ・ 全品揃っているか
- ・破損していないか

部品が見つからない場合や機器またはパッケージに破損が見られる場合は、直ちに運送業者に 連絡し、カスタマーサービスまでご連絡ください。 第3章 機器の詳細

3. 機器の詳細

3.1 説明

CELL LOCKER はポリカーボネート製チャンバーとメンブレンフィルター2枚で構成されています。 メンブレンフィルターは水蒸気と気体に対する透過性が高く、微生物に対して不透過性を示し ます。



☑ 1. CELL LOCKER

番号	名称
1	作業プレート
2	フィルタ付きメンブレン フレーム
3	ガスケット
4	ラッチ
5	トランスポートカバー (またはカバー)
6	ノブ



Makrolon 2528

図 2. 底のマーク

3.2 ご使用の前に

初めて使用する場合は、CELL LOCKER を清掃して殺菌してください。

留意 フィルター付きメンブレンフレームを挿入する前に、CELL LOCKER を清掃してください。

3.3 運転

CELL LOCKER は、6 ドア気密スクリーン付インキュベータ内に置いて、カバーを付けたままでも、 または外しても使用できます。

実験装置内に入れる際に、CELL LOCKERのトレイが開くか確認してください。



3.4 アプリケーションのバージョン

必要に応じて、消耗品キットを使用して CELL LOCKER のトランスポートカバーの付け外しが選択できます。



図 3. 消耗品キット

カバーは六角レンチを使用して作業トレイに取り付け可能で、移動する際の保護になります。



トランスポートカバー付き



トランスポートカバーなし



カバー付きのときの CELL LOCKER と培養容器の 取り出し



カバーなしのときの内扉の開放による培養 容器の取り出し



トランスポートカバーを摘まんでチャンバー 内のサンプルの出し入れ

図 4. CELL LOCKER のオプション

留意:

安全キャビネットや他のインキュ ベーターまたは研究室に移動する 場合は、保護のためにカバーを取 り付けてください。チャンバー内 でカバーを摘まんで開閉する際は、 両手を必要とする場合があります。 第4章 クリーニングと消毒

警告

4. クリーニングと消毒

作業開始の際に、以下の安全ガイドに従って実行してください。



健康被害の危険!



作業スペースの表面が汚染されていることがあります。汚染された洗浄液に接触す ると感染することがあります。消毒剤には有害物質が含まれていることがありま す。クリーニングや消毒を実行する際は、必ず安全上の注意と衛生管理のガイドラ インに従ってください。



セーフティゴーグルの着用!



保護手袋の着用!



呼吸用保護具の着用!

消毒剤の安全データシートや衛生管理者によって作成された安全上の注意に従って ください。

4.1 クリーニング

注意



非相容性洗浄剤!

CELL LOCKER はプラスチック製の部分があります。

溶剤を使用すると、プラスチックを侵す可能性があります。

強酸や強アルカリ溶液を使用すると、プラスチックの腐食を引き起こすことがあ ります。



注意 感湿性の部品!

> **CELL LOCKER**のメンブレンフィルターに洗浄剤をスプレーしないでください。メンブ レンフィルターのフレーム w を拭く際に、水分がメンブレンフィルターに接触しな いように注意してください。

ぬるま湯と中性洗剤の溶液で汚れや残留物、沈着物などを完全に除去します。

清潔な布と洗浄水で表面や作業トレイを拭きます。次に、清潔な布で表面の水気を拭き取りま す。 4.2 払拭消毒または噴霧消毒



アルコール消毒剤!

アルコール含有量が10%を超える消毒剤は、空気と混ざって可燃性や爆発性の高 い気体混合物を生成する可能性があります。

このような消毒剤を使用する場合は、消毒手順を行っている間、裸火や過度の熱 への露出を避けてください。

このような消毒剤を使用する場合は、必ず十分に換気された部屋で行ってください。 消毒剤で処理したら、消毒済の CELL LOCKER 部品を拭き、完全に水気を拭き取りま す。

アルコール含有消毒剤による火災・爆発の危険を避けるため、安全規定を順守し てください。



塩素系消毒剂!

注意

塩素系消毒剤は使用しないでください。

消毒剤はステンレス鋼に影響を与えないものを使用してください。



警告 除染または洗浄剤!

除染や洗浄剤の相容性についてご不明な点がありましたら、Thermo Electron LED GmbH か代理店までお問い合わせください。

- 1. すべてのサンプルを作業スペースから取り出し、安全な場所に保管します。
- 作業スペースと付属品の表面に消毒剤を噴霧するか、消毒剤で表面を拭き消毒します。メンブレンフィルターに消毒剤を噴霧しないようご注意ください。
- 3. 消毒剤の安全データシートに記載されている作用時間に従って放置します。
- 4. ぬるま湯に標準リンス剤を混ぜた液で、チャンバーの表面と作業トレイの消毒剤を拭き落 とします。温水で完全に洗い流します。
- 5. 高圧蒸気滅菌処理した水で洗浄した表面をさらに3~5回洗浄し、残留洗浄剤を完全に洗い 流します。
- 洗浄が完了したら、柔らかい滅菌布で浄化済の表面、エアバッフル、内部の水気を拭き取ります。

第4章 クリーニングと消毒

4.3 オートクレーブ

高速アスピレーターを使用した場合、CELL LOCKER は 121°C (250°F) で 15 分までのオート クレーブが可能です。高温でのオートクレーブ処理は、素材がそのために設計されていないた め、ダメージを受けるのでお勧めできません。



CELL LOCKER を高圧蒸気滅菌処理する前に、フィルター付きメンブレンフレームを取り外してください!

5. アフターサービス

保守作業はすべて記録しておきます。

5.1 メンブレンフィルターの交換

定期的にメンブレンフレームフィルターを確認し、少なくとも一年に一度交換してください。



1. 作業トレイを引いて取り出します。



2. 内側からメンブレンのフレームを押して取り外します。



3. メンブレンはビニール袋に入れ、高圧蒸気滅菌処理して有害廃棄物として処分します。



第5章 アフターサービス

4. 新しいメンブレンを取り出します。



- 5. 新しいメンブレンフレームを一方の半円部から挿入し、しっかりと押し込みます。
- ガスケットリップが正しくはめ込まれているかどうか目視で内部から確認してください。
 必要に応じて外側から押してください。



7.5と6に従って2つ目のメンブレンを挿入します。

8. 作業トレイを傾き止め上にはめ込み、内部に押し込みます。



第5章 アフターサービス

5.2 カバーガスケットの交換

カバーガスケットが完全な状態であるか、破損していないかなどについて定期的(例えば毎 月)に点検し、必要に応じて交換します。

1. 作業トレイを引いて取り出します。



2. カバーガスケットをノッチから取り外します。



ノッチに新しいガスケットをはめ込みます。ガスケットリップが外側になるように注意してください。





警告 汚染の危険!

破損したり、不適切に装着されたりしたメンブレンを使用すると、ユーザーや実験 室、環境などを汚染する恐れがあります。

4. 作業トレイを挿入し、内部に押し込みます。



6. 予備部品および摩耗部品

説明	部品番号	
メンブレンフレーム	50153148	
カバーシール (6 個)	50156433	
ステンレス製作業トレイ	50155808	
銅製作業トレイ	50155809	
消耗品キット	50161834	
Cell Locker カバー	50156434	

Thermo Fisher Scientific 承認済の原初スペアパーツと付属品のみが使用されている場合 (Thermo Fisher Scientific 非承認の第三者が製造したスペアが使用されている場合は、この 限定的保証は無効になります)。

7. 廃棄

警告 汚染の危険!

本装置は、感染性物質の作業や処理に使用されている場合がありますので、装置や 部品が汚染されている可能性があります。廃棄する前に、すべての装置の構成部品 を除染する必要があります。 ユニットの構成部品を完全に掃除してから、次に消毒・除染または高圧蒸気滅菌処 理を行ってください(使用したアプリケーションによる)。 処分する品目について、実施した除染処理の詳細が記載された異議な」宣言書を添

処分する品目について、実施した除染処理の詳細が記載された異議なし宣言書を添 付してください。

すべての装置の部品は、適切に汚染除去された後に廃棄することができます。 メンブレンフィルターは、必ず有害廃棄物の取り扱いに関する国内の法律および指令に基づい て処分してください。

構成部品の材質

構成部品	材質	
CELL LOCKER のハウジングおよびカバー	ポリカーボネート, Makrolon 2858	
ステンレス製作業トレイ	1.4301+2B / スチール 10088-2	
銅製作業トレイ	銅	
メンブレンフレーム	ABS 樹脂	
メンブレンフィルター	PALL社製 スーポア-200R PES メンブレン 0.2μm	

8. 仕様

説明	単位	数值
外形寸法(幅 × 高さ × 奥行)	mm	185 x 156 x 380
内部空間容積	I	約 5.4
	kg	1.8
作業トレイ1個当たりの積載量	kg	5
使用温度	°C/°F	4 - 55 / 39 - 131
高圧蒸気滅菌器サイクル		最大 12
高圧蒸気滅菌の温度	° C / ° F	最大 121 / 250



Find out more at thermofisher.com