thermo scientific



Thermo Scientific

STERI-CYCLE i160/ STERI-CYCLE i250 CO₂-Inkubator

Betriebsanleitung

50162762

März 2020

©2020 Thermo Fisher Scientific Inc. Alle Rechte vorbehalten.

Warenzeichen HERACELL VIOS™, Steri-run™, Steri-cycle™, iCan™, THRIVE™ und Cell locker™ sind eingetragene Warenzeichen von Thermo Scientific. Thermo Scientific ist eine Marke der Thermo Fisher Scientific Inc. Alle anderen in der Betriebsanleitung genannten Marken sind ausschließliches Eigentum der betreffenden Hersteller.

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

Die Thermo Electron LED GmbH ist einen Tochtergesellschaft von: Thermo Fisher Scientific Inc. 168 3rd Avenue Waltham, MA 02451 USA

Thermo Fisher Scientific Inc. stellt seinen Kunden dieses Dokument nach Erwerb eines Produktes für den Betrieb des Gerätes zur Verfügung. Dieses Dokument ist urheberrechtlich geschützt. Die Vervielfältigung - auch auszugsweise - ist ohne schriftliche Zustimmung von Thermo Fisher Scientific Inc. verboten.

Der Inhalt der Betriebsanleitung kann jederzeit und ohne Ankündigung geändert werden. Sämtliche in diesem Dokument enthaltenen technischen Angaben haben rein informatorischen Charakter und sind unverbindlich. Die in diesem Dokument enthaltenen Systemkonfigurationen und technischen Daten ersetzen etwaige frühere Angaben, die der Käufer erhalten hat.

Thermo Fisher Scientific Inc. erhebt keinen Anspruch auf die Vollständigkeit, Korrektheit und Fehlerfreiheit dieses Dokuments und haftet weder für hierin möglicherweise enthaltene Fehler oder Auslassungen noch für Folgeschäden, die sich aus der Verwendung dieses Dokuments ergeben, selbst wenn diese entsprechend den in diesem Dokument enthaltenen Angaben erfolgen sollte.

Dieses Dokument ist nicht Bestandteil eines Kaufvertrages zwischen Thermo Fisher Scientific Inc. und einem Käufer. Diese Dokument hat keinerlei Änderungseinfluss auf die Allgemeinen Verkaufsbedingungen, vielmehr haben die Allgemeinen Verkaufsbedingungen bei voneinander abweichenden Angaben in den Dokumenten in jedem Fall Vorrang.

Inhaltsverzeichnis

Vorwort		1
	Allgemeine Hinweise	1
	Kenndaten des Gerätes und der Gerätedokumentation	1
	Geräteidentifikation	1
	Einweisung des Bedienpersonals	1
	Gültigkeit der Anleitung	2
	Originaldokument für Übersetzung	2
	Gewährleistung	2
	Bedingungen Gewährleistung	2
	Erklärung der Sicherheitshinweise und Bildzeichen	3
	Sicherheitshinweise und Bildzeichen der Betriebsanleitung	3
	Ergänzende Bildzeichen zu den Sicherheitshinweisen:	4
	Zeichen am Gerät	5
	Verwendungszweck des Gerätes	5
	Bestimmungsgemäßer Gebrauch	5
	Bestimmungswidriger Gebrauch	6
	Normen und Richtlinien	6
	Sicherheitshinweise zu Gasen	7
	Einweisung des Personals:	7
	Sicherheitshinweise zu Kohlendioxid (CO2)	7
	Sicherheitshinweise zu Sauerstoff (O2)	8
	Sicherheitshinweise zu Stickstoff (N2)	8
Kapitel 1	Gerätelieferung	. 1-1
•	Verpackung	. 1-1
	Lieferkontrolle	. 1-1
	Lieferumfang Standardausstattung STERI-CYCLE i160 / i250	. 1-2
	Lieferumfang Zusatzausstattung STERI-CYCLE i160 / i250	. 1-2
	Lieferumfang Zusatzausstattung STERI-CYCLE i160	. 1-3
	Lieferumfang Zusatzausstattung STERI-CYCLE i250	. 1-3
Kapitel 2	Aufstellung und Installation	2-1
•	Umgebungsbedingungen	. 2-1
	Anforderungen:	. 2-1
	Raumbelüftung	. 2-2
	Platzbedarf	. 2-2
	Transport	. 2-3
	Geräte stapeln	. 2-3
	Stapelvarianten	. 2-7
	Untergestell mit Laufrollen montieren	. 2-8

Untergestell mit Laufrollen und Standfüßen (i160- i250-Ausführung)	
Nachrüst-/Umbauarbeiten	
Kapitel 3 Gerätebeschreibung	
STERI-CYCLE i160 / i250 Ansicht von vorne	
STERI-CYCLE i160 / i250 Rückansicht	
STERI-CYCLE i160 / STERI-CYCLE i250 mit durchgeschleifter Gasversorgung:	
Schutzeinrichtungen STERI-CYCLE i160 / i250	3-6
Nutzraum-Atmosphäre	
Temperatur:	
Relative Feuchte:	
Empfehlung zur Wasserqualität:	
Vorfilter	
HEPA-Filter und Luftführung	
CO2-Versorgung:	
O2-Zufuhr:	
N2-Zufuhr:	
Türschalter	
Sensorik	
Versorgungsschnittstelle	
Standardschnittstellen	
Optionale Schnittstellen	
Gasanschlüsse:	
Hinweisschild:	
USB-Schnittstelle:	
Schnittstelle 4-20 mA:	
Alarmkontakt:	
Netzanschluss:	
Komponenten des Nutzraumes	
Innenbehälter	
Glastür und optionale Gasblende	
Wasserreservoir	
Vorgehensweise beim Befüllen mit Wasser	
Heizsystem	
Rückwärtige Geräteöffnungen	
Regalsystem	
Elektromechanisches Türverschluss-Kit	
Kapitel 4 Inbetriebnahme	
• Gerät akklimatisieren	
Nutzraum aufbereiten	
Füllstandsanzeiger "MAX" und Vorfilter einsetzen	
Luftführung installieren	
HEPA-Filter und Wasserreservoirabdeckung einbauen	
Regalsystem installieren	
Tragprofile ein-/ausbauen	
Auflagebügel einsetzen	
Gerät nivellieren	
Geteilte Einlagebleche (optional) in STERI-CYCLE i250 einsetzen	4-11

	Gas anschließen	
	Gasdruckschläuche montieren	
	Gasanschluss ohne Gasmonitoring	
	CO2-Anschluss mit Gasmonitoring-System (optional)	
	Kombinierter CO2- und O2/N2-Anschluss mit Gasmonitoring-System (optional)	
	Netzanschluss	
	USB-Schnittstelle anschließen:	
	Alarmkontakt anschließen:	
	Schnittstelle 4-20mA anschließen:	
Kapitel 5	Betrieb	5-1
-	Gerät vorbereiten	
	Betrieb aufnehmen	
	Vorgehensweise beim Befüllen mit Wasser	
Kapitel 6	Bedienung	6-1
	Netzschalter	6-2
	Bedienfeld und Bedienstruktur	6-3
	Ausstattung ohne O2-/N2-Regelung:	6-4
	Ausstattung mit kombinierter CO2-/O2-/N2-Regelung (optional):	6-5
	Gliederung der Bedienebenen	6-5
	Werkseinstellungen der Regler des iCan™ Touchscreens	6-7
	Aufwärmphase der Sensoren der Regelkreise	6-7
	Verhalten der Tasten bei Einstellungen	6-8
	Temperatur-Sollwert einstellen	6-8
	CO2-Sollwert einstellen	6-9
	O2-Sollwert einstellen	6-10
	auto-start-Funktion	6-12
	auto-start aktivieren	6-14
	auto-start unterbrechen	6-15
	Steri-run aufrufen	6-16
	Benutzerkonfiguration	6-17
	Einstellungen/Setup	6-17
	Datenaufzeichnung	6-27
	Optionen	6-31
	Iconerklärung	6-40
	Tastensperre aktivieren/deaktivieren	6-42
	Softwareversionen	6-43
	Verlaufsanzeige skalieren	6-43
	Fehlermeldungen	6-45
	Reaktion auf das Ereignis Fehlermeldung	6-45
	Übertemperaturschutz zurücksetzen	6-46
	Maßnahmen nach einem Stromausfall	6-47
	Übersicht Fehlerursachen und Fehlerbeseitigung	6-48
Kapitel 7	Außerbetriebnahme	7-1
	Gerät außer Betrieb nehmen	7-1

Kapitel 8 Reinig	jung und Desinfektion	8-1
Rein	igung	
Dek	ontaminationsverfahren	
Desi	nfektion oder Steri-run vorbereiten	
Wise	ch-/Sprühdesinfektion	
Steri	-run-Dekontaminationsroutine	
A	blauf einer Steri-Run-Dekontaminationsroutine:	
St	eri-run aktivieren	
St	eri-run abbrechen	
St	eri-run Fehlerabbruch	
St	eri-run abschließen	
Ö	ffnen der Tür nach Abbruch der Steri-run-Dekontaminationsroutine	
Kapitel 9 Instan	dhaltung	
- Insp	ektionen und Kontrollen	
War	tungsintervalle	
Tem	peraturabgleich vorbereiten	
Tem	peraturabgleich durchführen	
CO2	2-Abgleich vorbereiten	
CO2	2-Abgleich durchführen	
HEI	PA-Filter tauschen	
Gase	einlassfilter tauschen	
Aust	ausch der Gerätesicherung	
Türe	lichtung wechseln	
Kapitel 10 Ents	orgung	
Übe	rsicht der verwendeten Materialien:	
Kanitel 11 Tech	nische Daten	11_1
STE	RI CVCI E :160	11.2
STE STE	RI-CYCLE 1160	
STE STE	RI-CYCLE 1100	
STE	RI-CYCLE 1250	
511		
Kapitel 12 Date	nkommunikation	
U	SB-Schnittstelle	
Т	reiber USB-Schnittstelle installieren	
Stru	ktur der Befehlssequenzen zur Datenkommunikation	
Be	eschreibung des Protokolls	
Übe	rsicht Allgemeine Parameter (Adressen 0xxx)	
A	llgemeine Parameter lesen	
Übe	rsicht Inkubator-Parameter (Adressen 2xxx)	
Pa	arameter (Basic) lesen	
Pa	arameter (interne Funktionen) lesen	
Stru	ktur Fehlerspeicher	
Sc	hema Datensatzstruktur Fehlerspeicher:	
Ü	bersicht der möglichen Fehlermeldungen in Hex-Codierung	
Ü	bersicht der möglichen Fehlermeldungen in Bit-Codierung	
	Allgemeiner Gerätestatus, Regelkreis Temperatur und CO2:	

Regelkreis O2 und Wasserstand:	
Struktur Datenlogger	
Schema Datensatzstruktur Datenlogger:	
Übersicht der möglichen Ereigniseinträge in Bit-Codierung	
Übersicht Ereigniseinträge Teil I:	
Übersicht Ereigniseinträge Teil II:	
Code-Beispiele Datenlogger	
Funktionen zur Abfrage Datenlogger	
Code-Beispiel zur Abfrage Datenlogger	
Programm STERI-CYCLE i160 / STERI-CYCLE i250	
STERI-CYCLE i160 & STERI-CYCLE i250 installieren	
STERI-CYCLE i160 & STERI-CYCLE i250 bedienen	
Kapitel 13 Gerätebuch	
Kapitel 14 Kontaktdaten Thermo Scientific	
Anhang A 6-fach-Gasblende für Cell Locker	

Anhang B Betriebsanleitung Cell Locker

Inhaltsverzeichnis

Abbildungsverzeichnis

Abbildung 2-1. Geräteabmessungen
Abbildung 2-2. Hebepunkte
Abbildung 2-3. Geräte stapeln2-4
Abbildung 2-4. Stapelrahmen und Untergestell mit Stapelelementen2-5
Abbildung 2-5. Verschrauben der Adapterplatte mit dem unteren Gerät2-5
Abbildung 2-6. Schrauben für die Trägerrahmen entfernen2-6
Abbildung 2-7. Trägerrahmen montieren2-6
Abbildung 2-8. Untergestell mit Laufrollen ohne Standfüße (i160-Ausführung)2-8
Abbildung 3-1. STERI-CYCLE i160 / STERI-CYCLE i250 Ansicht von vorne
Abbildung 3-2. STERI-CYCLE i160 / i250 Rückansicht
Abbildung 3-3. STERI-CYCLE i160 / STERI-CYCLE i250 mit durchgeschleifter Gasversorgung3-5
Abbildung 3-4. HEPA-Filter und Airbox
Abbildung 3-5. Luftführung
Abbildung 3-6. Türschalter
Abbildung 3-7. Sensoren für Temperatur, O2, CO2 und Luftfeuchte
Abbildung 3-8. Sensor für Wasserstand
Abbildung 3-9. Geräteschnittstellen (rechte Seite des Schaltkastens)
Abbildung 3-10. Optionale Geräteschnittstellen (linke Seite des Schaltkastens)
Abbildung 3-11. STERI-CYCLE i160 mit Glastür
Abbildung 3-12. STERI-CYCLE i160 mit 3-fach-Gasblende
Abbildung 3-13. Wasserreservoir
Abbildung 3-14. Befüll- und Ablassventil des Wasserreservoirs
Abbildung 3-15. Befüll- und Ablassventil des Inkubators
Abbildung 3-16. Befüllzylinder
Abbildung 3-17. Rückwärtige Geräteöffnungen
Abbildung 3-18. Komponenten des Regalsystems
Abbildung 3-19. Türverriegelungshebel und Notentriegelung an der Unterseite des Inkubators
Abbildung 4-1. Füllstandsanzeiger "MAX"
Abbildung 4-2. Füllstandsanzeiger "MAX" und Vorfilter
Abbildung 4-3. Luftführung zusammenbauen
Abbildung 4-4. Bauteile Luftführung STERI-CYCLE i160
Abbildung 4-5. HEPA-Filter und Airbox zusammenbauen
Abbildung 4-6. Airbox auf die Wasserreservoirabdeckung aufsetzen

A

Abbildung 4-7. Airbox einbauen	4-8
Abbildung 4-8. Ein-/Ausbau Regalsystem	4-9
Abbildung 4-9. Auflagebügel einsetzen	4-10
Abbildung 4-10. Einbau geteilter Einlagebleche	4-11
Abbildung 4-11. Gasdruckschläuche montieren	4-12
Abbildung 4-12. Gasanschluss ohne Gasmonitoring	
Abbildung 4-13. CO2-Anschluss mit Gasmonitoring-System (optional)	4-14
Abbildung 4-14. Kombinierter CO2- und O2-/N2-Anschluss mit Gasmonitoring-System (o	ptional)4-15
Abbildung 4-15. Netzanschluss	4-16
Abbildung 4-16. Beispiel Anschluss Alarmkontakt	4-19
Abbildung 4-17. Anschlussbelegung der Schnittstelle 4-20 mA	
Abbildung 5-1. Wasserreservoir	5-2
Abbildung 5-2. Befüll- und Ablassventil des Inkubators	5-3
Abbildung 5-3. Befüllzylinder	5-3
Abbildung 5-1. Füllstandsanzeiger "MAX"	5-4
Abbildung 6-1. Netzschalter	6-2
Abbildung 6-2. Hauptbildschirm: Tastsensitive Bildschirmbereiche	6-3
Abbildung 6-3. iCan [™] Touchscreen ohne Gasversorgung O2/N2	6-4
Abbildung 6-4. iCan™ Touchscreen mit kombiniertem Gasanschluss	6-5
Abbildung 6-5. Übersicht Menüstruktur	6-6
Abbildung 6-6. Anzeige Aufwärmphase	6-7
Abbildung 6-7. Temperatur-Anzeigefeld und Temperatur-Auswahlmenü	6-8
Abbildung 6-8. Temperatur-Sollwert einstellen	6-9
Abbildung 6-9. CO2-Anzeigefeld und CO2-Menü	6-9
Abbildung 6-10. CO2-Sollwert einstellen	6-10
Abbildung 6-11. O2-Anzeigefeld und O2-Menü	6-11
Abbildung 6-12. O2-Sollwert einstellen	6-11
Abbildung 6-13. auto-start aktivieren	6-14
Abbildung 6-14. Statusanzeige auto-start-Status	6-15
Abbildung 6-15. auto-start unterbrechen	6-16
Abbildung 6-16. Fehlermeldung nach Unterbrechen von auto-start	6-16
Abbildung 6-17. Menü Benutzerkonfiguration	6-17
Abbildung 6-18. Menü Einstellungen/Setup	6-18
Abbildung 6-19. Code für Tastensperre ändern	6-18
Abbildung 6-20. Code für Tastensperre ändern	6-19
Abbildung 6-21. Auswahlmenü Datum/Uhrzeit	6-20
Abbildung 6-22. Datum einstellen	6-20
Abbildung 6-23. Uhrzeit einstellen	6-21
Abbildung 6-24. Displayhelligkeit einstellen	6-22
Abbildung 6-25. Lautstärke des Tastentons einstellen	6-23

Abbildung 6-26. Auswahlmenü Schnittstelle USB	6-24
Abbildung 6-27. Baudrate USB-Schnittstelle einstellen	6-24
Abbildung 6-28. Sprache einstellen	6-25
Abbildung 6-29. Funktion für Erinnerungsintervall auswählen	6-26
Abbildung 6-30. Erinnerungsintervall für steri-run einstellen	6-27
Abbildung 6-31. Auswahlmenü Datenaufzeichnung	6-28
Abbildung 6-32. Ereignisse anzeigen	6-28
Abbildung 6-33. Speicherzyklus einstellen	6-29
Abbildung 6-34. Fehlertabelle einsehen	6-31
Abbildung 6-35. Auswahlmenü Optionen	6-32
Abbildung 6-36. Auswahlmenü Alarm	6-32
Abbildung 6-37. Alarmrelais einstellen	6-33
Abbildung 6-38. Low Feuchte einstellen	6-34
Abbildung 6-39. Gasblende einstellen	6-35
Abbildung 6-40. Wasserstandsensor einstellen	6-36
Abbildung 6-41. Alarmrelais einstellen	6-37
Abbildung 6-42. O2-Regelung ein-/ausschalten	6-38
Abbildung 6-43. HEPA-Konfiguration	6-39
Abbildung 6-44. HEPA-Filter aktivieren / deaktivieren	6-39
Abbildung 6-45. Iconerklärung	6-40
Abbildung 6-46. Icons Gasmonitoring	6-41
Abbildung 6-47. Tastensperre aktivieren/deaktivieren	6-43
Abbildung 6-48. Softwareversionen	6-43
Abbildung 6-49. Verlaufsanzeige für CO2-Konzentration aufrufen	6-44
Abbildung 6-50. Verlaufsanzeige für CO2-Konzentration aufrufen	6-44
Abbildung 6-51. Ereignis Fehlermeldung	6-46
Abbildung 6-52. Fehlermeldung Übertemperatur	6-46
Abbildung 6-53. Fehlermeldung Übertemperatur	6-47
Abbildung 7-1. Befüll- und Ablassventil des Wasserreservoirs	7-1
Abbildung 8-1. Befüll- und Ablassventil des Wasserreservoirs	8-2
Abbildung 8-2. HEPA-Filter und Airbox	8-5
Abbildung 8-3. Luftführung	8-6
Abbildung 8-4. Phasen der Dekontaminationsroutine	8-9
Abbildung 8-5. Menü steri-run - Ablauf	8-9
Abbildung 8-6. Steri-run aktivieren	8-10
Abbildung 8-7. Steri-run unterbrechen, abbrechen	8-11
Abbildung 8-8. Steri-run beenden	8-12
Abbildung 8-9. Türverriegelungshebel und Notentriegelung an der Unterseite des Inkubators	8-13
Abbildung 9-1. Temperaturabgleich vorbereiten	9-3
Abbildung 9-2. Temperatur-Anzeigefeld und Temperatur-Auswahlmenü	9-4

Abbildung 9-3. Temperaturabgleich durchführen	9-4
Abbildung 9-4. Messöffnung Gasblenden	9-6
Abbildung 9-5. CO 2-Abgleich durchführen	9-7
Abbildung 9-6. Airbox ausbauen	9-8
Abbildung 9-7. HEPA-Filter einbauen	9-9
Abbildung 9-8. Gaseinlassfilter montieren	9-10

Vorwort

Allgemeine Hinweise

Kenndaten des Gerätes und der Gerätedokumentation Geräteidentifikation

Geräteidentifikation

Gerätebezeichnung: Gerätebezeichnung: CO₂-Inkubator *STERI-CYCLE i160 STERI-CYCLE i250*

Zuordnung der Produktdokumentation Dokumentations-CD 50143971

Zertifizierung und Qualitätsaudit:

Konformität:CE-KonformitätskennzeichnungPrüfzeichen:TÜV GS, cCSAus, EAC

Einweisung des Bedienpersonals

Diese Betriebsanleitung beschreibt den CO_2 -Inkubator **STERI-CYCLE i160 / STERI-CYCLE i250**. Der CO_2 -Inkubator ist nach dem aktuellen Stand der Technik gefertigt und wurde vor der Auslieferung auf einwandfreie Funktion geprüft. Dennoch können von diesem Gerät Gefahren ausgehen. Vor allem dann, wenn es von nicht ausreichend unterwiesenem Personal bedient wird oder wenn es unsachgemäß, und nicht wie für den bestimmungsgemäßen Gebrauch vorgesehen, verwendet wird. Zur Unfallverhütung sollten deshalb die folgenden Verfahrensweisen berücksichtigt werden:

- Der CO₂-Inkubator darf nur von eingewiesenem und autorisiertem Personal bedient werden.
- Für Personal, das mit diesem Gerät arbeitet, sind auf der Grundlage dieser Betriebsanleitung, der gültigen Sicherheitsdatenblätter, der betrieblichen Hygienerichtlinien und der entsprechenden Technischen Regeln vom Betreiber schriftliche Verfahrensanweisungen zu erstellen, insbesondere:
 - welche Dekontaminationsmaßnahmen für das Gerät und die verwendeten Hilfsmittel anzuwenden sind,
 - welche Sicherheitsmaßnahmen bei der Verwendung von Gasen und Druckgasbehältern einzuhalten sind,
 - welche Maßnahmen bei Unfällen zu ergreifen sind.
- Instandsetzungsarbeiten am Gerät dürfen nur von ausgebildetem und autorisiertem Fachpersonal ausgeführt werden.

Gültigkeit der Anleitung

Originaldokument für Übersetzung

- Der Inhalt dieser Betriebsanleitung kann jederzeit und ohne Ankündigung geändert werden.
- Für Übersetzungen in Fremdsprachen ist die englische Fassung dieser Betriebsanleitung verbindlich.
- Bewahren Sie die Betriebsanleitung in der Nähe des Gerätes auf, damit jederzeit Sicherheitshinweise und wichtige Informationen zur Bedienung nachgeschlagen werden können.

Bei Fragen, die nach Ihrer Auffassung in dieser Betriebsanleitung nicht ausführlich genug behandelt werden, wenden Sie sich zu Ihrer eigenen Sicherheit an Thermo Scientific.

Gewährleistung

Bedingungen Gewährleistung

Thermo Scientific gewährleistet die Sicherheit und die Funktionstüchtigkeit des CO2-Inkubators (für 2 Jahre) nur unter der Bedingung, dass:

- das Gerät ausschließlich für den bestimmungsgemäßen Gebrauch eingesetzt wird und gemäß den Angaben dieser Betriebsanleitung bedient und instand gehalten wird,
- keine baulichen Veränderungen am Gerät vorgenommen werden,
- nur originale und von Thermo Scientific zugelassene Ersatzteile oder Zubehör verwendet werden,
- Inspektionen und Wartungsarbeiten entsprechend den vorgegebenen Zeitintervallen durchgeführt werden.

Der Gewährleistungszeitraum der Cell Locker und Zubehörteile beträgt 1 Jahr. Die Gewährleistungszeitraum beginnt mit der Auslieferung des Gerätes an den Besteller.

Erklärung der Sicherheitshinweise und Bildzeichen

Sicherheitshinweise und Bildzeichen der Betriebsanleitung



WARNUNG Bezeichnet eine Gefahrensituation, die zu Tod oder schwerer Verletzung führen könnte, wenn sie nicht vermieden wird.



VORSICHT Bezeichnet eine Gefahrensituation, die zu leichter oder mäßiger Verletzung führen könnte, wenn sie nicht vermieden wird.



VORSICHT Bezeichnet eine Situation, die zu Sachschäden führen kann, wenn sie nicht vermieden wird.

Hinweis Gibt Anwendungstipps und nützliche Informationen.

Ergänzende Bildzeichen zu den Sicherheitshinweisen:



Schutzhandschuhe tragen!



Schutzbrille tragen!



Gefährliche Flüssigkeiten!



Stromschlag!



Heiße Oberflächen!



Feuergefahr!



Explosionsgefahr!



Erstickungsgefahr!

Zeichen am Gerät



CE-Konformitätskennzeichnung: bestätigt die Konformität gemäß den Richtlinien der EU



TÜV-geprüfte Sicherheit



Prüfzeichen USA/Kanada



EAC-Konformitätskennzeichnung: bestätigt die Konformität mit sämtlichen technischen Vorschriften der Eurasischen Zollunion (Russland, Kasachstan und Weißrussland)



Betriebsanleitung beachten!



Heiße Oberflächen!

Verwendungszweck des Gerätes

Bestimmungsgemäßer Gebrauch

Dieser CO_2 -Inkubator ist für die Aufbereitung und Kultivierung von Zellkulturen konzipiert, ist jedoch nicht für medizinische Anwendungen oder für die In-vitro-Diagnose zu verwenden. Zu diesem Zweck werden im Arbeitsraum des Gerätes durch exakte Regelung folgender Parameter kontrollierte physiologischen Umgebungsbedingungen hergestellt:

- Temperatur
- CO₂-Gehalt
- O₂/N₂-Gehalt
- Relative Feuchte

Typischerweise wurde der $\rm CO_2$ -Inkubator für Installation und Betrieb in folgenden Anwendungsbereichen entwickelt:

- Labore für zellbiologische und biotechnologische Versuche der Sicherheitsstufen L1, L2 und L3
- Medizinisch-mikrobiologische Labore gemäß DIN EN 12128
- Forschungslabore in Kliniken und Krankenhäusern

Dieses Gerät ist ausschließlich für den professionellen Einsatz konzipiert und darf nur von entsprechend geschultem Fachpersonal betrieben werden.

Bestimmungswidriger Gebrauch

Es dürfen keine Zell- und Gewebekulturen im Gerät verwendet werden, die nicht den Bestimmungen der Sicherheitsstufen L1, L2 und L3 entsprechen. Es dürfen keine Gewebe, Stoffe oder Flüssigkeiten als Proben verwendet werden, die:

- die leicht entflammbar oder explosiv sind,
- deren freigesetzte Dämpfe in Verbindung mit Luft brennbare oder explosive Gemische bilden,
- die Gifte freisetzen.

Der Cell Locker dürfen nicht in Medizingeräten verwendet werden.

Normen und Richtlinien

Das Gerät entspricht den Sicherheitsanforderungen folgender Normen und Richtlinien:

- Niederspannungsrichtlinie 2014/35/EG
- IEC 61010-1:2010+Korr. 2011, Sicherheitsbestimmungen für elektrische Mess-, Steuer-, Regelund Laborgeräte, Teil 1: Allgemeine Anforderungen
- IEC 61010-2-010:2003, Sicherheitsbestimmungen für elektrische Mess-, Steuer-, Regel- und Laborgeräte, Teil 2-010: Besondere Anforderungen an Laborgeräte für das Erhitzen von Stoffen
- EMV-Richtlinie 2014/30/EU
- IEC 61326-1:2012, Elektrische Mess-, Steuer-, Regel- und Laborgeräte EMV-Anforderungen, Teil 1: Allgemeine Anforderungen
- FCC 47, Part 15, § 15.107, EMV-Regulierung FCC
- FCC 47. Part 15, § 15.109, EMV-Regulierung FCC

In anderen Ländern sind die entsprechenden nationalen Vorschriften bindend.

Sicherheitshinweise zu Gasen

Hinweis Installationsarbeiten:

Arbeiten an Zuleitungen und Druckgasbehältern, Flaschen oder Sammelbehältern, in denen CO_2 bzw. O_2/N_2 bereitgehalten wird, dürfen nur von Fachpersonal mit den dazu geeigneten Werkzeugen durchgeführt werden.

Hinweis

Es ist sicherzustellen, dass die Arbeitsplatzgrenzwerte für CO $_2$ bzw. O $_2/\rm N_2$ nicht überschritten werden.

Für die Bundesrepublik Deutschland wird empfohlen die TRGS 900 zu beachten, in anderen Ländern können abweichende Grenzwerte gelten. Die länderspezifischen, nationalen Arbeitsplatzgrenzwerte sind einzuhalten.

Einweisung des Personals:

Personal, das an Geräten mit CO_2 -Zufuhr arbeitet, ist vor Aufnahme der Tätigkeit über den besonderen Umgang mit CO_2 zu unterweisen:

- die sachgemäße Bedienung von Druckgasbehältern und Gasversorgungsanlagen,
- die Meldepflicht von Schäden und Mängeln an den CO₂-Zuleitungen,
- Maßnahmen, die bei Unfällen und Störungen zu treffen sind.

Die Unterweisungen sind in angemessenen Zeitabständen zu wiederholen. Die besonderen Bedienungsanweisungen des Gaslieferanten sind in die Unterweisung einzubeziehen.

Sicherheitshinweise zu Kohlendioxid (CO₂)

 CO_2 wird als gesundheitsgefährdendes Gas eingestuft. Deshalb müssen bei Inbetriebnahme und während der Nutzung des CO_2 -Inkubators folgende Sicherheitsmaßnahmen eingehalten werden:



Erstickungsgefahr!

Wird Kohlendioxid (CO₂) in großen Mengen in die Raumatmosphäre freigesetzt, besteht Erstickungsgefahr.

Bei Austritt von CO2 sofort die Sicherheitsmaßnahmen einleiten!

- Sofort den Raum verlassen und die Zugänge sichern!
- Sicherheitsdienst oder Feuerwehr benachrichtigen!

Sicherheitshinweise zu Sauerstoff (O₂)

O2 ist ein verbrennungsförderndes Gas und reagiert explosiv mit fetthaltigen Materialien.

Sauerstoffexplosion!



Sauerstoff (O_2) reagiert explosiv mit Ölen, Fetten und Schmiermitteln. Bei Kontakt des hochverdichteten Sauerstoffs mit fett- und ölhaltigen Substanzen besteht Explosionsgefahr! Bedien- und Anzeigefolie!

• Zum Reinigen dieser Geräteteile nur öl- und fettfreie Reinigungsmittel benutzen.

Alle Anschlüsse und Komponenten der Sauerstoffanlage frei von öl-, fett- und schmiermittelhaltigen Stoffen halten!

Feuergefahr!



Austretender Sauerstoff (O_2) ist extrem verbrennungsfördernd. Kein offenes Feuer in der Umgebung von sauerstoffführenden Anlagen benutzen!

• In der Nähe von Sauerstoffanlagen nicht rauchen.

Die Komponenten der Sauerstoffanlage keiner starken Hitzeeinwirkung aussetzen.

Sicherheitshinweise zu Stickstoff (N₂)

Stickstoff mischt sich leicht mit Luft. Hohe Konzentrationen von Stickstoff verringern den Sauerstoffgehalt in der Luft.



Erstickungsgefahr!

Wird Stickstoff (N_2) in großen Mengen in die Raumatmosphäre freigesetzt, besteht Erstickungsgefahr wegen Sauerstoffmangels. Bei Austritt von N_2 die Sicherheitsmaßnahmen einleiten!

- Sofort den Raum verlassen und die Zugänge sichern!
- Sicherheitsdienst oder Feuerwehr benachrichtigen!

Gerätelieferung

Inhalt

- "Verpackung" auf Seite 1-1
- "Lieferkontrolle" auf Seite 1-1
- "Lieferumfang Standardausstattung STERI-CYCLE i160 / i250" auf Seite 1-2
- "Lieferumfang Zusatzausstattung STERI-CYCLE i160 / i250" auf Seite 1-2
- "Lieferumfang Zusatzausstattung STERI-CYCLE i250" auf Seite 1-3

Verpackung

Der CO₂-Inkubator **STERI-CYCLE i160** / **STERI-CYCLE i250** wird in einer stabilen Verpackungskiste geliefert. Sämtliche Verpackungsmaterialien können getrennt werden und sind wiederverwertbar:

- Verpackungskarton: Altpapier
- Kunststoffschaumteil: Styropor (FCKW-frei)
- Verpackungsfolie: Polyethylen
- Verpackungsbänder: Polypropylen
- Füße: Polypropylen
- Palette: unbehandeltes Holz

Lieferkontrolle

Überprüfen Sie sofort nach Anlieferung des Gerätes:

- die Vollständigkeit der Lieferung,
- den Lieferzustand des Gerätes.

Ist die Lieferung unvollständig oder sind Transportschäden an Gerät und Verpackung, insbesondere Schäden durch Feuchtigkeit und Wasser, feststellbar, informieren Sie umgehend den Spediteur, sowie den technischen Support.

Lieferumfang Standardausstattung STERI-CYCLE i160 / i250

Mitgelieferte Gerätekomponenten	Stückzahl
Wasserreservoirabdeckung	1
Max. Wasserstandsensor	1
Luftleitblech Rückwand	1
Luftleitblech Decke	1
Airbox incl. Dichtung	1
Vorfilter	1
Einlagebleche	3
Tragprofile für Einlagebleche	4
Auflagebügel für Einlagebleche	6
Stopfen für Rohrdurchführung	1
Netzkabel	1
CO ₂ -Anschlussschlauchsatz	1
Betriebsanleitung	1
Schnellverschluss mit Schlauch für Wasserablauf	1

Lieferumfang Zusatzausstattung STERI-CYCLE i160 / i250

Lieferumfang Zusatzausstattung Sauerstoffregelung	Stückzahl
O ₂ -Sensorkopf	1
O ₂ -Anschlussschlauchsatz	1
Lieferumfang Zusatzausstattung Gasflaschenmonitor CO ₂	
Anschlussschlauchsatz CO ₂ -Gasflaschenmonitor	1
Lieferumfang Zusatzausstattung Gasflaschenmonitor O ₂ /N ₂	
Anschlussschlauchsatz O ₂ /N ₂₋ Gasflaschenmonitor	1
Lieferumfang Zusatzausstattung Filter	
HEPA-Filter oder VOC-Filter	1
Lieferumfang Zusatzausstattung Schnittstelle 420 mA	
Stecker 420mA	1

Lieferumfang Zusatzausstattung STERI-CYCLE i160

Komponenten	Stückzahl
3-fach-Gasblende	1
6-fach-Gasblende	1
Cell Locker	6
Befüllzylinder	1

Lieferumfang Zusatzausstattung STERI-CYCLE i250

Lieferumfang 250 Liter Geräte mit geteilten Einlagen	Stückzahl
Trägerrahmen	3
Geteilte Einlagebleche	6
6-fach-Gasblende	1

1 Gerätelieferung Lieferumfang Zusatzausstattung STERI-CYCLE i250

2

Aufstellung und Installation

Inhalt

- "Umgebungsbedingungen" auf Seite 2-1
- "Raumbelüftung" auf Seite 2-2
- "Platzbedarf" auf Seite 2-2
- "Transport" auf Seite 2-3
- "Geräte stapeln" auf Seite 2-3
- "Nachrüst-/Umbauarbeiten" auf Seite 2-11

Umgebungsbedingungen

Das Gerät darf nur an Aufstellungsorten betrieben werden, welche die unten aufgeführten besonderen Umgebungsbedingungen erfüllen:

Anforderungen:

- Zugluftfreier und trockener Aufstellungsort.
- Die Mindestabstände zu angrenzenden Flächen nach allen Seiten müssen eingehalten werden, siehe "Platzbedarf" auf Seite 2-2.
- Der Betriebsraum muss mit einer geeigneten Raumlüftung ausgestattet sein.
- Feste, ebene und nicht brennbare Aufstellfläche.
- Verwenden Sie dazu einen tragfähigen, vibrationsfreien Unterbau (Untergestell, Labortisch), welcher der Belastung durch die Gerätegewichte und Zuladung standhält (besonders beim Stapeln der Geräte).
- Das Gerät ist für einen Betrieb an einem Standort auf max. 2000 m über NN ausgelegt.
- Um eine gängige, konstante Inkubationstemperatur von 37° C zu erreichen, muss die Raumtemperatur im Bereich von +18° C bis +34° C liegen.
- Relative Luftfeuchte bis max. 80 %.
- Keine direkte Sonneneinstrahlung.
- Es dürfen keine Geräte mit hoher Wärmeabstrahlung in naher Umgebung des **STERI-CYCLE i160** / **STERI-CYCLE i250** installiert oder abgestellt sein.

Raumbelüftung

Beim Einleiten von CO_2 - $/O_2$ - $/N_2$ entsteht ein geringfügiger Überdruck im Nutzraum des Inkubators, der durch die Druckausgleichsöffnung in den Betriebsraum abströmt und so ausgeglichen wird. Durch den Druckausgleich und durch das Öffnen der Glastür/Gasblende im laufenden Betrieb wird in sehr geringen Mengen CO_2 - $/O_2$ - $/N_2$ in den Betriebsraum abgegeben. Die Raumlüftung muss das ausströmende Gas gefahrlos ins Freie abführen können.

Zudem kann im Dauerbetrieb durch die Energieabgabe des Gerätes eine Veränderung des Raumklimas entstehen.

- STERI-CYCLE i160 / STERI-CYCLE i250 deshalb nur in ausreichend belüfteten Räumen aufstellen.
- Das Gerät nicht in unbelüfteten Nischen aufstellen.
- Die Raumlüftung sollte eine technische Lüftung sein, die den Vorgaben der nationalen Richtlinien für Laboratorien entspricht, oder eine entsprechend leistungsstarke Lüftungsanlage.

Platzbedarf



WARNUNG NOT-AUS!

Die Steckdose für den Netzanschluss muss für den Fall eines NOT-AUS immer zugänglich sein. Der Netzstecker muss für den Benutzer klar erkennbar und jederzeit frei zugänglich sein.



Abbildung 2-1. Geräteabmessungen

Hinweis Zugänglichkeit der Geräte:

Für eine bessere Zugänglichkeit bei Pflege- und Wartungsarbeiten wird empfohlen, größere seitliche und größere rückwärtige Abstände einzuhalten. Zum Ausschwenken der Tür muss bei Geräten mit 6-fach-Gasblende und Cell Locker an der Türanschlagseite ein Abstand von >= 31 cm / 12,2" zur Wand vorhanden sein.

Transport

Zum Transport darf das Gerät nicht an den Türen oder Anbauteilen, wie z. B. dem rückwärtigen Schaltkasten, angehoben werden.



Abbildung 2-2. Hebepunkte

Hinweis Hebepunkte:

Das Gerät nur an den in der Abbildung gekennzeichneten Hebepunkten belasten.

VORSICHT Schwere Lasten! Vorsicht beim Anheben!



Zur Vermeidung überbelastungsbedingter Verletzungen, wie z.B. Zerrungen und Bandscheibenschäden, niemals versuchen, den Inkubator allein anzuheben! Zur Vermeidung von Verletzungen durch herabfallende Lasten sind beim Anheben des Inkubators stets persönliche Schutzausrüstungen zu tragen, wie z.B. Sicherheitsschuhe. Zur Vermeidung von Quetschungen der Finger oder Hände (insbesondere durch Einklemmen beim

Schließen der Tür) oder Beschädigung des Inkubators dürfen ausschließlich die in der vorstehenden Abbildung dargestellten Hebepunkte benutzt werden.

Geräte stapeln

STERI-CYCLE i160 / **STERI-CYCLE i250** ist zum Stapeln von maximal 2 Geräten desselben Gerätetyps geeignet. Dazu wird zwischen beiden Geräten ein optionaler Stapeladapter (1/Abbildung 2-3) eingesetzt.

Als Option ist ein bewegliches Untergestell (2/Abbildung 2-3) zum Verfahren der Geräte lieferbar. Die Rollen des Untergestelles sind mit einer Feststellbremse ausgestattet.

Hinweis

Bei der Installation der Stapeladapterplatte und der Stapelung der Geräte ist die dem Stapeladapter beiliegende Montageanleitung zu beachten.

Hinweis Verfahren von gestapelten Geräten:

Vor dem Verfahren von gestapelten Geräten ist sicherzustellen, dass die Trägerrahmen entfernt wurden!

Gestapelte Geräte auf einem beweglichen Untergestell sind ausschließlich in Räumen mit glattem Fußboden ohne Gefälle zu verschieben.



Abbildung 2-3. Geräte stapeln

1. Unteres Gerät mit den Gerätefüßen (3/Abbildung 2-4) auf die Stapelelemente (1/Abbildung 2-4 an der Oberseite des beweglichen Untergestells (2/Abbildung 2-4) aufsetzen.



Abbildung 2-4. Stapelrahmen und Untergestell mit Stapelelementen

- 2. Adapterplatte (7/Abbildung 2-5) mit der Unterseite auf die Gerätedecke des unteren Gerätes (8/Abbildung 2-5) setzen.
- 3. Die Bohrungen in der Anschlusslasche (9/Abbildung 2-5) der Adapterplatte (7/Abbildung 2-5) auf beiden Seiten mit den Bohrungen in der Rückwand des unteren Gerätes (8/Abbildung 2-5) fluchten.





- 4. Adapterplatte (7/Abbildung 2-5) mit den mitgelieferten Befestigungsschrauben mit der Rückwand des unteren Gerätes (8/Abbildung 2-5) verschrauben.
- 5. Adapterplatte (7/Abbildung 2-5) an der anderen Ecke des Geräts ebenfalls mit zwei Schrauben befestigen.
- 6. Das zu stapelnde Gerät mit den Gerätefüßen (6/Abbildung 2-5) auf die Stapelelemente (1/Abbildung 2-4) an der Oberseite der Adapterplatte aufsetzen.

Die Geräte werden durch das Eigengewicht der Adapterplatte (7/Abbildung 2-5) und des oberen Gerätes (10/Abbildung 2-5) aufeinander fixiert.

7. Die vier Schrauben an der oberen hinteren Kante des oberen Gerätes entfernen, siehe Pfeile in Abbildung 2-6.



Abbildung 2-6. Schrauben für die Trägerrahmen entfernen

- 8. Beide Enden der Trägerrahmen (4/Abbildung 2-7) um einen Winkel von ca. 90 ° verbiegen, um zwei rechtwinklige Laschen zu erhalten.
- 9. Die vier Schrauben verwenden, um den Trägerrahmen an der Geräterückseite zu befestigen, wobei die Lasche nach unten auf die Geräteseite zeigt.
- 10. Sicherstellen, dass die Stapelfüße des oberen Gerätes korrekt mit dem Stapeladapter ausgerichtet sind.
- 11. Die Trägerrahmen an einem tragfähigen Gebäudeteil mit geeignetem Montagematerial, wie Holzschrauben und Dübel, befestigen.



Abbildung 2-7. Trägerrahmen montieren



VORSICHT Gefahr des Kippens und Herunterfallens gestapelter Geräte! Um die Gefahr des Kippens des Gerätes zu vermeiden, muss der Trägerrahmen an einer Gebäudewand montiert werden, die der Last der gestapelten Geräte standhalten kann. Die Installation darf nur durch qualifiziertes Personal durchgeführt werden. Für die Befestigung der Trägerrahmen am Gebäudeteil geeignete Schrauben und Dübel mit einer Tragfähigkeit von mindestens 25 kg verwenden.



VORSICHT Transport gestapelter Geräte!

Die Stapelelemente sind keine Verbindungselemente. Gestapelte Geräte auf einem beweglichen Untergestell sind ausschließlich in Räumen mit glattem Fußboden ohne Gefälle zu verschieben.

Hinweis Fixierung der beweglichen Untergestelle:

Werden die Geräte auf beweglichen Untergestellen aufgestellt, muss sichergestellt sein, dass die Rollen während des Betriebes der Inkubatoren mit einer Feststellbremse fixiert sind und dass die Rollen wegen der höheren Standsicherheit nach vorne ausgerichtet sind.

Kondensatbildung bei Betrieb mit gestapelten Geräten:

Generell muss beim Betrieb von gestapelten Geräten vom Typ **STERI-CYCLE i160** und **i250** eine Adapterplatte, als thermische Trennung, verwendet werden.

Werden gestapelte Geräte bei einer Umgebungstemperatur von mehr als 28 °C betrieben, tritt, während das untere Gerät die Steri-run-Dekontaminationsroutine durchläuft, am oberen Gerät ein Übertemperaturfehler auf. Dann kann es am oberen Gerät zu Kondensatbildung kommen.

Stapelvarianten

Mögliche Stapelkombinationen		Untere Stapelposition				
		HERAcell Vios 160i	HERAcell Vios 250i	HERAcell 150i	HERAcell 240i	
		Steri-Cycle i160	Steri-Cycle i250			
lposition	HERAcell Vios 160i	50148171	50154522	50148172		
	Steri-Cycle i160					
òtape	HERAcell Vios 250i		50154522		50148175	
Obere (Steri-Cycle i250					

Weitere Informationen finden Sie in der Montageanweisung des Stapeladapter-Satzes.

Untergestell mit Laufrollen montieren





Abbildung 2-8. Untergestell mit Laufrollen ohne Standfüße (i160-Ausführung)



VORSICHT Gewährleistung der Stabilität Sicherstellen, dass die Laufrollen des Untergestells während des Gerätebetriebs verriegelt sind und dass die Laufrollen nach vorne zeigen.

- Das Untergestell mit dem Hebel [3] an der Oberseite der Laufrollen verriegeln.
- Die Gerätefüße [1] exakt auf die 4 Stapelplättchen[2] platzieren. Zum Positionieren darf das Gerät nicht an den Türen oder Anbauteilen, wie z. B. dem rückwärtigen Schaltkasten, angehoben werden.

Untergestell mit Laufrollen und Standfüßen (i160- i250-Ausführung)

Lieferumfang

1 Untergestell, 6 Schrauben M4x16



• Die 6 Schrauben unten an der Rückseite des Gerätes entfernen. Sie werden durch die sechs neuen Schrauben ersetzt.

- Das Gerät auf das Untergestell platzieren und dabei die korrekte Positionierung der Standfüße achten. Sicherstellen, dass die Bohrungen der entfernten Schrauben exakt mit den Bohrungen in den Ecken des Untergestells ausgerichtet sind.
- Die 6 neuen Schrauben einführen und festschrauben.



• Das Untergestell und das aufgesetzte Gerät vorsichtig zum gewünschten Aufstellungsort schieben.

Hinweis

Die Standfüße mit einem Schraubenschlüssel festschrauben, bevor das Gerät am Untergestell verfahren wird.

• Das Untergestell durch Festschrauben der Standfüße mit einem Schraubenschlüssel einstellen, bis dieses waagerecht steht.



VORSICHT Gewährleistung der Stabilität

Sicherstellen, dass die Standfüße herausgeschraubt und mit einem Schraubenschlüssel korrekt eingestellt wurden, nachdem das Gerät am gewünschten Aufstellungsort positioniert wurde.



Hinweis

Das Gerät nur an den in der Abbildung gekennzeichneten Hebepunkten anheben.



VORSICHT Schwere Lasten! Vorsicht beim Anheben!

Zur Vermeidung überbelastungsbedingter Verletzungen, wie z.B. Zerrungen und Bandscheibenschäden, niemals versuchen, den Inkubator allein anzuheben! Zur Vermeidung von Verletzungen durch herabfallende Lasten sind beim Anheben des Inkubators stets persönliche Schutzausrüstungen zu tragen, wie z. B. Sicherheitsschuhe. Zur Vermeidung von Quetschungen der Finger oder Hände (insbesondere durch Einklemmen beim Schließen der Tür) oder Beschädigung des Inkubators dürfen ausschließlich die in der vorstehenden Abbildung dargestellten Hebepunkte benutzt werden.



VORSICHT

Kippen des Gerätes auf einem Untergestell mit Laufrollen.



Beim Verfahren des Gerätes ist größte Vorsicht geboten. Durch unvorsichtiges Abbremsen oder Beschleunigen des Gerätes kann dieses vom Untergestell mit Laufrollen kippen. Das Gerät nicht mit geöffneter Fronttür verfahren. Das Untergestell mit Laufrollen sollte nur in der Nähe des Aufstellungsortes im Labor verwendet werden und nicht zu Transportzwecken.

Teilenummer	Beschreibung
50145394	Untergestell mit Laufrollen für HERAcell VIOS 160i und Steri-Cycle i160
50154551	Untergestell mit Laufrollen für HERAcell VIOS 160i und Steri-Cycle i160
50154407	Untergestell mit Laufrollen für HERAcell VIOS 250i und Steri-Cycle i250

Nachrüst-/Umbauarbeiten

Eine Nachrüstung der Standardversion ist möglich für folgende Bauteile:

STERI-CYCLE i160

- 3-fach-Gasblende (die Gasblende wird anstelle der Glastür montiert),
- Türanschlagwechsel Außentür und Glastür,
- 6-fach-Gasblende mit Cell Locker,
- abschließbare Außentür,
- Infrarot (IR) CO₂-Sensor (ersetzt den serienmäßigen WLD-Sensor),
- Messdatenschnittstelle 4-20 mA.
- Gasflaschenmonitor

STERI-CYCLE i250

- 3-fach-Gasblende (die Gasblende wird anstelle der Glastür montiert),
- Türanschlagwechsel Außentür und Glastür,
- abschließbare Außentür,
- geteilte Einlagen,
- Infrarot (IR) CO₂-Sensor (ersetzt den serienmäßigen WLD-Sensor),
- Messdatenschnittstelle 4-20 mA.
- Gasflaschenmonitor

Hinweis Umbauarbeiten:

Nachrüst- und Umbauarbeiten dürfen nur vom Technischen Service der Thermo Electron LED GmbH durchgeführt werden. **2** Aufstellung und Installation Nachrüst-/Umbauarbeiten
Gerätebeschreibung

Inhalt

- "STERI-CYCLE i160 / i250 Ansicht von vorne" auf Seite 3-2
- "STERI-CYCLE i160 / i250 Rückansicht" auf Seite 3-4
- "Schutzeinrichtungen STERI-CYCLE i160 / i250" auf Seite 3-6
- "Nutzraum-Atmosphäre" auf Seite 3-6
- "Türschalter" auf Seite 3-10
- "Sensorik" auf Seite 3-11
- "Versorgungsschnittstelle" auf Seite 3-13
- "Komponenten des Nutzraumes" auf Seite 3-17

STERI-CYCLE i160 / i250 Ansicht von vorne



Abbildung 3-1. STERI-CYCLE i160 / STERI-CYCLE i250 Ansicht von vorne

- 1. Druckausgleichsöffnung /Rohrdurchführung mit Stopfen
- 2. Türschalter für Glastür
- 3. Temperatursensor
- 4. O₂-Sensor (optional)
- 5. Türgriff Glastür bzw. Gasblende
- 6. Glastür bzw. Gasblende (optional)
- 7. Gerätetür
- 8. Magnettürdichtung, austauschbar
- 9. Vorfilter
- 10. Wasserablauf
- 11. Fuß, höhenverstellbar
- 12. Dichtung, Glastür, austauschbar
- 13. Airbox mit Dichtung und HEPA-Filter

14. -

- 15. Tragprofile
- 16. Einlageblech mit Auflagebügel
- 17. -
- 18. IR CO₂-Sensor
- 19. Türgriff Außentür
- 20. iCan[™] Touchscreen (Bedienteil)
- 21. Netzschalter
- 22. a = Elektromechanischer Türverschluss / b Verschlusshaken an der Außentür (nur an Geräten mit optionalem Türverschluss)

STERI-CYCLE i160 / i250 Rückansicht



Abbildung 3-2. STERI-CYCLE i160 / i250 Rückansicht

- 1. Schaltkasten mit Versorgungsschnittstellen für kombinierten Gasanschluss (optional) CO_2 und O_2/N_2 ohne Gasmonitoring
- 2. Schema: Gasanschluss CO2 und O2/N2 ohne Gasmonitoring
- 3. O₂/N₂-Gasflasche
- 4. CO₂-Gasflasche
- 5. Primäre CO₂-Gasversorgung A bei optionalem Gasmonitoring
- 6. Sekundäre CO₂-Gasversorgung B bei optionalem Gasmonitoring
- 7. Schaltkasten mit Versorgungsschnittstellen für kombinierten Gasanschluss (optional) CO_2 und O_2/N_2 mit Gasmonitoring (optional)
- 8. Primäre O₂-/N₂-Gasversorgung A bei optionalem Gasmonitoring
- 9. Sekundäre O₂-/N₂-Gasversorgung B bei optionalem Gasmonitoring
- 10. Schema: Gasanschluss CO_2 und O_2/N_2 mit optionalem Gasmonitoring

11. Verteileranschluss zur Weiterschleifung der CO₂-Gasversorgung für 3 weitere Geräte (unabhängig vom Gerätetyp)

Verdeckt: Verteileranschluss zur Weiterschleifung der O₂-/N₂-Gasversorgung für 3 weitere Geräte, auf gegenüberliegender Seite des Schaltkastens

STERI-CYCLE i160 / STERI-CYCLE i250 mit durchgeschleifter Gasversorgung:



Abbildung 3-3. STERI-CYCLE i160 / STERI-CYCLE i250 mit durchgeschleifter Gasversorgung

- 1. Schaltkasten mit Versorgungsschnittstellen für kombinierten Gasanschluss (optional) CO_2 und O_2/N_2 mit Gasmonitoring (optional)
- 2. Verteileranschluss zur Weiterschleifung der CO₂-Gasversorgung für 3 weitere Geräte (unabhängig vom Gerätetyp)
- 3. Verteileranschluss (verdeckt) zur Weiterschleifung der O₂-/N₂-Gasversorgung für 3 weitere Geräte (unabhängig vom Gerätetyp)
- 4. T-Stück zum Verbinden der Druckschläuche
- 5. Druckschlauch zum Weiterschleifen der Gasversorgung

- 6. Schema: Gasanschluss CO₂ und O₂/N₂ mit Gasmonitoring (optional) für CO₂ und O₂
- 7. **STERI-CYCLE i160** / **STERI-CYCLE i250** mit durchgeschleifter Gasversorgung; dazu ist nur ein kombinierter Gasanschluss (optional) CO_2 und O_2/N_2 erforderlich.

Schutzeinrichtungen STERI-CYCLE i160 / i250

Das Gerät ist mit folgenden Schutzeinrichtungen ausgerüstet:

- Ein Türschalter unterbricht beim Öffnen der Glastür die Nutzraumheizung und die Gaszufuhr für CO₂/O₂/N₂.
- Bei Einsatz der Option Gasmonitoring schaltet ein Gaswächter die Gasversorgung auf eine gefüllte Gasflasche um.
- Ein Übertemperaturschutz schützt die Proben im Fehlerfall vor einer schädlichen Überhitzung.
- Eine Druckausgleichsöffnung sorgt für den Druckausgleich im Nutzraum.
- Die Schaltung des Alarmrelais, akustische und optische Warnsignale weisen auf Fehler während des Betriebes hin.

Nutzraum-Atmosphäre

Im Nutzraum des Inkubators werden die besonderen physiologischen Umgebungsbedingungen für die Aufbereitung und Kultivierung von Zell- und Gewebekulturen simuliert. Die Nutzraum-Atmosphäre wird dabei von folgenden Faktoren bestimmt:

- Temperatur
- Relative Feuchte
- CO₂-Konzentration
- O₂-Konzentration (optional)

Temperatur:

Für einen störungsfreien Betrieb muss die Temperatur des Betriebsraumes mindestens 18 °C betragen und die Inkubationstemperatur mindestens 3 °C über der Raumtemperatur des Betriebsraumes liegen. Das Heizsystem regelt die Inkubationstemperatur von dieser Temperaturschwelle bis 55 °C. Die Beheizung des Innenbehälters mit unabhängigen Heizkreisen und die zusätzliche separate Beheizung der Außentür sorgen dafür, dass sich an den Seitenwänden und der Decke des Nutzraumes sowie an der Glastür bzw. Gasblende kein Kondensat bildet.

Relative Feuchte:

Die Beheizung des Nutzraumes fördert die Verdunstung des Wassers und sorgt so für eine konstante Feuchte im Nutzraum. Für den laufenden Betrieb muss eine ausreichende Menge an aufbereitetem Wasser mit der empfohlenen Wasserqualität vorrätig gehalten werden:

• Maximale Füllmenge für STERI-CYCLE i160 und STERI-CYCLE i250: 3 l.

Empfehlung zur Wasserqualität:

Um einen störungsfreien Betrieb zu gewährleisten, muss das Wasserreservoir mit sterilisiertem und destilliertem oder entsprechend aufbereitetem Wasser befüllt werden. Die Leitfähigkeit des Wassers sollte im Bereich von 1 bis 20 µS/cm liegen (der elektrische Widerstand im Bereich von 50 kOhmcm bis 1 MOhmcm).



VORSICHT Erlöschen der Gewährleistung!

Bei der Verwendung von chloriertem Leitungswasser oder Wasserzusätzen, die Chlor enthalten erlischt die Gewährleistung. Die Gewährleistung erlischt ebenso bei der Verwendung von Reinstwasser (Ultrapure Water), dessen Leitfähigkeit außerhalb des Bereiches von 1 bis 20 μ S/cm liegt und dessen elektrischer Widerstand außerhalb des Bereiches 50 kOhmcm bis 1 MOhmcm liegt. Bei Fragen bitte den Technischen Service von Thermo Fisher Scientific kontaktieren.

VORSICHT Kein Trink- oder Reinstwasser im Feuchtigkeitsbehälter



Für die Verwendung im integrierten Wasserreservoir wird steriles destilliertes Wasser oder qualitativ entsprechend behandeltes Wasser empfohlen. Die zulässige Leitfähigkeit sollte im Bereich von 1 bis 20 μS/cm liegen (der elektrische Widerstand sollte im Bereich von 50 K-Ohm-cm bis 1 M-Ohm-cm liegen). Der pH-Wert sollte zwischen 7 und 9 betragen. Reinstwasser Typ 1 oder entionisiertes (DI) Wasser mit einem elektrischen Widerstand von annähernd oder gleich 18,2 M-Ohm-cm enthält sehr wenig Ionen und zieht daher aktiv Ionen aus den innen liegenden Komponenten, wodurch es zu Beschädigungen von Edelstahl, Kupfer und Glas kommen kann. Ist ausschließlich DI- oder Typ 1-Wasser erhältlich, kann eine sterile Lösung aus schwachem Natriumbikarbonat hinzugefügt werden, um den pH-Wert zu erhöhen und Ionen hinzuzufügen (empfohlen wird: 84 mg/l (1 mmol/l).

VORSICHT Keine chloridhaltigen Desinfektionsmittel verwenden!

Obwohl Edelstahl korrosionsfest ist, ist er nicht korrosionsbeständig. Viele Chemikalien haben einen negativen Effekt auf Edelstahl, insbesondere Chlor und Derivate mit oxidierender Wirkung. Die Zugabe von chloridhaltigen Desinfektionsmitteln oder Kupfersulfat in das Wasser als konstantes Desinfektionsmittel wird nicht empfohlen, da diese das verbindende Abflussgelenk, welches aus einer Stahl/Kupfer-Legierung besteht, beschädigen können.Zur Reinigung des Innenraums wird zum Ausspülen eine milde Wasser-Seifenlösung empfohlen, um Rückstände zu entfernen. Wischen Sie die Innenflächen und -teile mit einem verdünnten quatären Ammonium-Desinfektionsmittel ab. Anschließend mit 70 % Alkohol abwischen, um alle verbleibenden Spuren des Desinfektionsmittels zu entfernen.

Unter normalen Betriebsbedingungen stellt sich bei einer üblichen Inkubationstemperatur von 37 $^{\circ}$ C im Nutzraum eine konstante, relative Feuchte von ca. 93 % ein.

Bildet sich aufgrund der hohen relativen Feuchte an den Kulturgefäßen ein Tauniederschlag, kann die Feuchte im Nutzraum auf einen niedrigeren Wert angepasst werden. Mit der Aktivierung der Low Feuchte wird die relative Feuchte im Nutzraum von ca. 93 % auf ca. 90 % abgesenkt. Die Veränderung beansprucht eine lange Anpassungsphase. Damit sie wirksam eine Betauung von Kulturgefäßen verhindert, muss sie als Dauereinstellung angewendet werden.

Eine Anleitung zum Aktivieren der Funktion Low Feuchte ist im Abschnitt "Low Feuchte einstellen" auf Seite 6-33 enthalten.

Vorfilter

Im vorderen Bereich der Wasserreservoirabdeckung ist ein Vorfilter aufgesetzt. . Das Vorfilter besteht aus einem doppelten Metalldrahtgeflecht mit Silikonrahmen und ist autoklavierbar sowie temperaturbeständig. Während der Steri-run-Dekontaminationsroutine muss das Vorfilter im Gerät eingebaut bleiben, zum Befüllen des Wasserreservoirs wird es herausgenommen.

HEPA-Filter und Luftführung

Der Luftstrom vom Wasserreservoir in Richtung Nutzraum wird zur Minimierung des Kontaminierungsrisikos durch ein HEPA-Filter geleitet. Das Filter arbeitet mit einem Abscheidegrad von 99,998 % bezogen auf eine Partikelgröße von 0,3 µm (HEPA-Filterqualität).

Das HEPA-Filter (2/Abbildung 3-4) wird von unten in die Airbox (1/Abbildung 3-4) eingesetzt. Die Airbox sitzt auf einem Sockel auf der Wasserreservoirabdeckung (2/Abbildung 4-6) und wird an den Lüftereinlass geschoben.



Abbildung 3-4. HEPA-Filter und Airbox

Eine Anleitung zum Aktivieren der Überwachung für den HEPA-Filter ist im Abschnitt "HEPA-Filter aktivieren / deaktivieren:" auf Seite 6-39 enthalten.

Die Luftführung kanalisiert den Luftstrom vom Lüfter entlang der Rückwand (3/Abbildung 3-5) zur Decke des Nutzraums und sorgt für eine optimale Temperaturverteilung. Gleichzeitig leitet sie die einströmenden Prozessgase in den Nutzraum ein und sorgt für eine optimale Durchmischung des Gasgemischs.



Abbildung 3-5. Luftführung

Die Luftführung besteht aus zwei Blechprofilen:

- 1. Deckenkanal (1/Abbildung 3-5)
- 2. Rückwandkanal (2/Abbildung 3-5)
- 3. Rückwand des Nutzraums (3/Abbildung 3-5)

Luftführung und HEPA-Filter lassen sich ohne Werkzeug ein- und ausbauen.

CO₂-Versorgung:

Um die Wachstumsbedingungen der Zell- und Gewebekulturen sicherzustellen, wird dem Nutzraum CO₂ zugeführt.

Der pH-Wert in den Bikarbonat-gepufferten Kulturmedien wird wesentlich vom CO₂-Gehalt der Nutzraumatmosphäre beeinflusst.

Der CO₂-Gehalt der Nutzraumatmosphäre kann im Bereich von 0 - 20 % geregelt werden. Das zugeführte CO₂ muss eines der folgenden Qualitätsmerkmale aufweisen:

- Reinheit min. 99,5 %
- medizinische Gasqualität.

O₂-Zufuhr:

Soll der CO₂-Inkubator mit mehr als 21 % Sauerstoff betrieben werden, wird dem Nutzraum Sauerstoff zugeführt (optional).

Der O2-Gehalt der Nutzraumatmosphäre kann im Bereich von 21 % - 90 % geregelt werden.

Bei Betrieb mit höheren Sauerstoffkonzentrationen sind die Brandschutzhinweise im Vorwort im Abschnitt Sicherheitshinweise zu Gasen auf Seite 7 zu beachten.

N₂-Zufuhr:

Soll der Sauerstoffgehalt im Betrieb unter 21 % (Luftsauerstoffgehalt) gesenkt werden, wird dem Nutzraum Stickstoff zugeführt. Der O₂-Gehalt der Nutzraumatmosphäre kann so, abhängig von der Ausführung des Sensors, geregelt werden.

Türschalter



Abbildung 3-6. Türschalter

An der Oberkante der Nutzraumöffnung ist ein Türschalter installiert. Wird der Türschalter durch Öffnen der Glastür aktiviert, werden die Gaszufuhr und die Beheizung des Nutzraumes unterbrochen. Am Bedienfeld wird eine Hinweismeldung angezeigt.

Ist die Tür länger als 30 s geöffnet, ertönt ein kurzer Signalton. Bleibt die Tür länger als 10 min geöffnet, wird ein Hupton ausgelöst und das Alarmrelais geschaltet.

Die Außentür kann nur verschlossen werden, wenn die Glastür korrekt verriegelt ist.

Hinweis Ausführung mit Gasblende:

Bei Geräten mit optionaler Gasblende wird die oben beschriebene Funktion des Türschalters bereits beim Öffnen der Außentür aktiviert.

Sensorik

In der Rückwand des Nutzraums sind das Lüfterrad und die Sensormodule installiert:

- Sensor zur Erfassung der Nutzraumtemperatur und des Übertemperaturschutzes (1/Abbildung 3-7).
- O₂-Sensor (optional) zur Erfassung des Sauerstoffgehaltes der Nutzraumatmosphäre (2/Abbildung 3-7).
- CO₂-Sensor zur Erfassung des CO₂-Gehaltes der Nutzraumatmosphäre (3/Abbildung 3-7). Je nach Ausstattung ist an dieser Position ein WLD-Sensor oder IR-Sensor (Option) eingebaut.
- rH-Sensor zur Erfassung der Luftfeuchte der Nutzraumatmosphäre (4/Abbildung 3-7). Der rH-Sensor (Option) dient der Feuchtekompensation in Verbindung mit dem WLD-Sensor und ist im Betrieb mit IR-Sensor nicht vorhanden.



Abbildung 3-7. Sensoren für Temperatur, O2, CO2 und Luftfeuchte

 Der Wasserstand-Sensor (1/Abbildung 3-8) alarmiert den Anwender, wenn das Wasserreservoir (2/Abbildung 3-8) nachgefüllt werden muss. Wenn das Wasserreservoir bis auf 0,5 l aufgebraucht ist, wird in der Displayanzeige im Feld rH die Meldung Fehler - rH - Kein Wasser (siehe auch "Fehlermeldungen" auf Seite 6-45) eingeblendet.



Abbildung 3-8. Sensor für Wasserstand

• Zusätzlich ist eine mechanische/optische Anzeige für den maximalen Wasserstand als Hilfe zum Befüllen vorhanden (siehe/Abbildung 3-8).

Der Sensor zur Messung der Nutzraumtemperatur sowie der CO₂-Sensor und der optionale O₂-Sensor sind Teil des Regelsystems des Gerätes. Die von ihnen gelieferten Messwerte werden mit den eingestellten Sollwerten verglichen. Das Regelsystem regelt auf dieser Datengrundlage die Heizung und die Gaszufuhr für CO₂-/O₂-/N₂.

Der Lüfter sorgt für eine Durchmischung der eingeleiteten Gase mit der angefeuchteten Luft und somit für eine gleichmäßige Temperaturverteilung im Nutzraum.

Der Übertemperaturschutz ist werkseitig vorprogrammiert und nur durch geschultes Service-Personal veränderbar. Er schützt die eingelagerten Kulturen vor Überhitzung. Wird die Soll-Temperatur um mehr als 1 °C überschritten, wird der Übertemperaturschutz aktiviert und die Nutzraumtemperatur automatisch auf den eingestellten Sollwert abgesenkt. Der Inkubationsbetrieb wird somit auch im Fehlerfall fortgesetzt. Jede Aktivierung des Übertemperaturschutzes löst gleichzeitig auch ein optisches Warnsignal aus. Ist der Übertemperaturschutz aktiviert, werden:

- eine Fehlermeldung (Temp.-Istwert über) und ein Hupton ausgegeben,
- das Alarmrelais geschaltet.

Wird die Fehlermeldung quittiert, wird als Hinweis auf die Aktivierung des Übertemperaturschutzes das Icon Übertemperatur am Display angezeigt und das Temperatur-Anzeigefeld ist rot hinterlegt.

Versorgungsschnittstelle

Standardschnittstellen



Abbildung 3-9. Geräteschnittstellen (rechte Seite des Schaltkastens)

Sämtliche Versorgungsanschlüsse sind an der Versorgungsschnittstelle (Schaltkasten) an der Geräterückseite installiert.

Auf der rechten Seite (Abbildung 3-9) des Schaltkastens sitzen die Anschlüsse der Grundausstattung sowie einige optionale Elemente:

- 1. Vier Anschlüsse an den internen Gasverteiler zum Betrieb mit der Option Gasmonitoring sowie zum Durchschleifen der CO₂-Gasversorgung an weitere Geräte
- 2. O₂-Anschlussstutzen (nicht vorhanden bei CO₂- und O₂-/N₂-Gasversorgung mit Option Gasmonitoring)
- 3. CO₂-Anschlussstutzen
- 4. Hinweisschild
- 5. USB-Schnittstelle
- 6. Alarmkontakt
- 7. Netzanschluss

Optionale Schnittstellen



Abbildung 3-10. Optionale Geräteschnittstellen (linke Seite des Schaltkastens)

Die Anschlüsse auf der linken Seite (Abbildung 3-10) sind nur bei Geräten mit entsprechenden Optionen vorhanden:

- 1. Vier Anschlüsse an den internen Gasverteiler zum Betrieb mit der Option Gasmonitoring sowie zum Durchschleifen der ebenfalls optionalen O₂-Gasversorgung an weitere Geräte
- 2. Schnittstelle 4...20 mA (Option)
- 3. O₂-Anschlussstutzen (Option; nur vorhanden bei CO₂- und O₂-/N₂-Gasversorgung mit Option Gasmonitoring)

Gasanschlüsse:

Die Gaszuleitungen zwischen Gerät und Gasversorgungsanlage werden über die mitgelieferten Anschlussschläuche hergestellt. CO_2 und O_2/N_2 werden dem Gerät über Anschlussstutzen (2 und 3/Abbildung 3-9 sowie 3/Abbildung 3-10) zugeführt.

Alle Prozessgase müssen am Gerät mit einem fest voreingestellten, nicht veränderbarem Druck im Bereich von min. 0,8 und max. 1,0 bar zugeführt werden.

Vor dem Einleiten in den Nutzraum durchströmen die Gase ein Gaseinlassfilter mit einem Abscheidegrad von 99,998 % bezogen auf eine Partikelgröße von 0,3 µm (HEPA-Filterqualität). Die Abbildung zeigt den kombinierten Gasanschluss (optional) ohne das Gasmonitoring-System (optional).

Hinweisschild:

Das Hinweisschild (4/Abbildung 3-9) enthält Angaben zur Versorgung mit Gasen, zur Belegung der Alarmkontaktanschlüsse und zur elektrischen Absicherung des Gerätes.

USB-Schnittstelle:

Über die USB-Schnittstelle (5/Abbildung 3-9) kann der Inkubator mit einem PC verbunden werden. Diese Verbindung - USB 1.1 / USB 2.0 / USB 3.0 full speed compatible - ermöglicht den schnellen (auch temporären) Zugang zu den wichtigsten Betriebsparametern (Temperatur, CO_2 -/ O_2 -/ N_2 -Konzentration, Fehlercodes, etc.).

Schnittstelle 4-20 mA:

Die optionale Schnittstelle 4-20 mA (2/Abbildung 3-10) wandelt die angezeigten digitalen Istwerte der Temperatur, CO_2 -Konzentration und O_2 -Konzentration (optional) in einen kontinuierlichen Ausgangsstrom von jeweils 4-20 mA um. Hierzu wird ein Mikroprozessor in Verbindung mit einem 4-Kanal-D/A-Wandler eingesetzt. Der D/A-Wandler besitzt eine Auflösung von 16 Bit pro Kanal, entsprechend 65536 Schritten.

An die Schnittstelle können kundenseitig verschiedene externe Messgeräte mit 4-20 mA-Eingängen angeschlossen werden, z.B. das Thermo Scientific Smart-Vue Wireless Monitoring System (gehört nicht zum Lieferumfang).

Kanal Nr.	Ausgegebener Parameter (Istwert)	Auflösung	Messbereich	Ausgangs- signal
1	Temperatur	0,08 MA/°C	0200 °C	420mA
2	CO ₂ -Konzentration	0,8 mA/%	020% CO ₂	420mA
3	O ₂ - Konzentration*	0,16 mA/%	0100% O ₂	420mA
4	frei			

Die Ausgangssignale sind in der nachfolgenden Tabelle aufgeführt:

Die mit * gekennzeichneten Parameter werden nur ausgegeben, wenn die entsprechende Optionen vorhanden ist.

Die Messwerte folgen der Funktion:

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

 $I_0 = 4m A$, MW = Messwert (°C, % oder rH%) $I_{max} = 20m A$, MB = Messbereich

Die Bedeutung der Ausgangssignale wird in der nachfolgenden Tabelle erklärt:

Strom	Bedeutung
4 mA	Der Messwert ist kleiner oder gleich dem Minimalwert des Messbereiches.
20 mA	Der Messwert ist größer oder gleich dem Maximalwert des Messbereiches.
2 mA	Das Gerät ist voll funktionsfähig, aber es liegt kein gültiger Wert vor (z.B. der optionale Sensor ist nicht vorhanden, das Gerät befindet sich in der Anwärmphase, o.ä.).
0 mA	Fehler

Die vier 4-20 mA-Ausgangssignale werden auf einen steckbaren, 8-poligen Anschlussblock (3,5 mm) herausgeführt. Der entsprechende Gegenstecker gehört zum Lieferumfang. Die Pinbelegung kann der nachfolgenden Tabelle entnommen werden.

Pin Nr.	Kanalbezeichnung
1	Kanal 1: Temperatur
2	Masse Kanal 1
3	Kanal 2: CO2
4	Masse Kanal 2
5	Kanal 3: O2
6	Masse Kanal 3
7	Kanal 4: frei
8	Masse Kanal 4

Alarmkontakt:

Das Gerät kann an ein kundenseitiges, externes Meldesystem (z.B. Telefonanlage, Gebäudeleitsystem, optische oder akustische Alarmmelder) angeschlossen werden.

Dazu ist ein potentialfreier Alarmkontakt im Gerät vorinstalliert. Dieser Kontakt wird am Schaltkasten auf der Geräterückseite herausgeführt (6/Abbildung 3-9).

Hinweis Alarmkontakt:

Der Alarmkontakt schaltet bei allen von den Regelkreisen gemeldeten Fehlern (siehe Abschnitt Fehlermeldungen).

Netzanschluss:

Der Netzanschluss des Gerätes erfolgt mit einem Kabel mit Kaltgerätestecker an der Kaltgerätesteckdose im Schaltkasten (7/Abbildung 3-9). Der Netzstecker muss für den Benutzer klar erkennbar und jederzeit frei zugänglich sein.

Komponenten des Nutzraumes

Der Nutzraum des Inkubators ist so konzipiert, dass durch technische Maßnahmen zur Vermeidung von Kondenswasserbildung und ein direkt im Nutzraum eingebautes HEPA-Filter-System keine störende Kontaminierung auftreten kann. Das HEPA-Filter-System im Nutzraum schützt das zur Luftbefeuchtung verwendete Wasser, sorgt für eine reinraumähnliche Luftqualität (Reinraumklasse ISO 5) und ist dabei völlig frei von Rückwirkungen auf den Lagerbereich der Kulturen.

Innenbehälter

Der Nutzraum des Inkubators ist so ausgelegt, dass alle den Inkubationsbetrieb beeinträchtigenden Kontaminationen ausgeschlossen werden. Dies wird durch Unterbinden der Kondensatbildung und den Einsatz eines in den Nutzraum eingebauten HEPA-Filtrierungssystems erreicht, das ohne Abstriche bei der zum Inkubieren der Kulturen nutzbaren Fläche den zum Befeuchten verwendeten Wasservorrat schützt und für eine Reinraum-Luftqualität nach ISO 5 sorgt.

Materialien des Innenbehälters

Die Standardversion ist ausgestattet mit einem Innenbehälter aus Edelstahl oder Kupfer.

Je nach Werkstoff des Innenbehälters sind Nutzraumkomponenten wie Luftführung und Regalsystem ebenfalls aus demselben Edelstahl- oder Kupfermaterial gefertigt.

Die Airbox für den HEPA-Filter ist aus temperaturbeständigem Kunststoff gefertigt und muss auch eingesetzt bleiben, wenn die Steri-run-Dekontaminationsroutine durchlaufen wird.



WARNUNG Der HEPA-Filtereinsatz ist nur bis 60 °C temperaturbeständig, nicht autoklavierbar und muss vor der Steri-run-Dekontamination entfernt werden.

Hinweis Oxidation der Kupferteile:

Durch die Einwirkung von Hitze und Luftfeuchtigkeit oxidiert das Kupfermaterial des Innenbehälters. Die Kupferteile verfärben sich deshalb bereits während des Testbetriebes bei der Geräteprüfung.

Die Oxidationsschicht sollte bei der routinemäßigen Reinigung nicht entfernt werden, weil auf ihr die antimikrobielle Wirkung des Kupfermaterials beruht.

Die Komponenten der Einlegeböden, die Airbox, die Luftführung und die Abdeckung des Wasserreservoirs lassen sich einfach ohne Werkzeugeinsatz herausnehmen, so dass zur Reinigung und manuellen Desinfektion des Gerätes lediglich ein einfach zu behandelnder, flächenreduzierter Innenbehälter bleibt.

Glastür und optionale Gasblende

In der Standardversion sind die Geräte **STERI-CYCLE i160** und **STERI-CYCLE i250** mit einer einfach verriegelten Tür aus Einscheibensicherheitsglas ausgerüstet.



Abbildung 3-11. STERI-CYCLE i160 mit Glastür

Geräte, die mit der optionalen Gasblende ausgerüstet sind, haben ein vielfach geringeres Kontaminationsrisiko wegen der wesentlich kleineren Öffnungsquerschnitte beim Zugriff auf die Proben und erreichen kürzere Erholungszeiten der Inkubationsparameter für:

- Nutzraumtemperatur,
- CO₂-Konzentration
- O₂-/N₂-Konzentration,
- relative Feuchte

Gasblenden sind als optionale Ausstattungsvarianten für beide Gerätetypen verfügbar:

- STERI-CYCLE i160: 3-fach-Gasblende (siehe Abbildung 3-12),
- STERI-CYCLE i250: 6-fach-Gasblende.
- Optionale 6-fach-Gasblende mit Cell Locker (siehe Anhang).



Abbildung 3-12. STERI-CYCLE i160 mit 3-fach-Gasblende

Jede Glastür der Gasblende (2/Abbildung 3-12) besitzt eine eigene Verriegelung (1/Abbildung 3-12).

Wasserreservoir

Das Wasserreservoir ist in den Boden des Innenbehälters integriert und durch eine Abdeckung vom Nutzraum getrennt. Ein Wasserablass (2/Abbildung 3-13) im vorderen Bereich des Wasserreservoirs ermöglicht eine schnelle Entleerung über das Befüll- und Ablassventil (3/Abbildung 3-13) an der Gerätevorderseite.

Im vorderen Bereich der Wasserreservoirabdeckung ist ein Vorfilter (4/Abbildung 3-13) aufgesetzt. Das Vorfilter besteht aus einem doppelten Metalldrahtgeflecht mit Silikonrahmen und ist autoklavierbar sowie temperaturbeständig. Während der Steri-run-Dekontaminationsroutine muss das Vorfilter im Gerät eingebaut bleiben, zum Befüllen des Wasserreservoirs wird es herausgenommen. Während der Steri-run-Dekontaminationsroutine muss das Vorfilter im Gerät eingebaut bleiben, zum Befüllen des Wasserreservoirs wird es herausgenommen.

Weitere Informationen zum Befüllen mit der 6-fach-Gasblende sind im Anhang zu finden.



Abbildung 3-13. Wasserreservoir

Das Wasserreservoir wird durch den im Abschnitt "Sensorik" auf Seite 3-11 beschriebenen Wasserstandsensor überwacht.

Als Markierung für die maximale Füllmenge ist ein Füllstandsanzeiger mit der Beschriftung "MAX" (1/Abbildung 3-13) in das Wasserreservoir eingehängt. Als maximale Füllmenge für das Wasserreservoir sind 3 l zulässig.

Um Störungen der Nutzraumatmosphäre beim Wassertausch im laufenden Inkubationsbetrieb so gering wie möglich zu halten, verfügt das Gerät über einen frontseitigen Schnellablass. Durch Einsetzen des mitgelieferten Ablassschlauchs in die frontseitige Schnellablassarmatur wird die sofortige Entleerung eingeleitet.



Abbildung 3-14. Befüll- und Ablassventil des Wasserreservoirs

Vorgehensweise beim Befüllen mit Wasser



Abbildung 3-15. Befüll- und Ablassventil des Inkubators

Der Inkubator kann mit einem Befüllzylinder über das Wasserventil erneut mit Wasser befüllt werden.



Abbildung 3-16. Befüllzylinder

Die Haken der Befüllvorrichtung in die Verschlüsse der Glastüren einhängen (siehe oben). Den Schlauch an das Befüll- und Ablassventil anschließen.

Der Befüllzylinder bestimmt nach dem Prinzip der kommunizierenden Röhren die Höhe des Wasserstands im Inkubator. Die minimale Füllmenge beträgt 0,5 l, die maximale Füllmenge beträgt 3,5 l. Der Zylinder enthält ca. 0,4 l Wasser.

- Den Zylinder in den Wasserablass einsetzen. Einige Zeit warten, bis sich der Wasserstand stabilisiert hat, sodass dieser abgelesen werden kann.
- Um das Befüllen durch die Einwirkung der Schwerkraft zu ermöglichen, Zylinder bis zum Maximum mit sterilem destilliertem Wasser befüllen. Dieser Vorgang muss ggf. mehrmals wiederholt werden.
- Wasser hinzufügen, bis die maximale Füllstandsanzeige erreicht ist.

Heizsystem

Zur Beheizung des Nutzraumes dient ein Flächenheizsystem. Die Heizelemente sind so angeordnet, dass eine Kondensatbildung oberhalb des Wasserreservoirs weitestgehend verhindert wird. Die Außentür des Gerätes und der Umfang der Türöffnung wird ebenfalls beheizt. Die Abstrahlung der Wärme auf die innenliegende Glastür/Gasblende verhindert einen Kondensatniederschlag. Die Sicht in den Nutzraum des Gerätes bleibt trotz hoher Feuchte immer frei.

Rückwärtige Geräteöffnungen



Abbildung 3-17. Rückwärtige Geräteöffnungen

Die durch einen Stopfen (1/Abbildung 3-17)verschließbare Rohrdurchführung mit Einsatz (Ø Abbildung 3-1742 mm) ermöglicht die Verlegung von Leitungen, Schläuchen oder zusätzlichen Sensoren in den Nutzraum des Gerätes.

Die Druckausgleichsöffnung (3/Abbildung 3-17) unterhalb der Rohrdurchführung an der Rückwand des Gerätes sorgt für den Druckausgleich zwischen Gerätenutzraum und Betriebsraum.

Hinweis

Um Schäden am Silikon zu vermeiden, entfernen Sie unbedingt den Silikonstopfen aus dem Innenbehälter und stecken Sie ihn in die Außenseite der Zugangsöffnung, bevor Sie mit der Hochtemperatur-Dekontamination beginnen.

Hinweis Betriebsbedingungen:

Beim Betrieb von Hilfsgeräten im Nutzraum des CO_2 -Inkubators müssen die Anforderungen an die Umgebungsbedingungen beachtet werden (vgl. Tabelle). Die in den Nutzraum eingebrachte Energie hat Einfluss auf den Beginn des Regelbereiches der Temperatur. Beim Einbringen zusätzlicher Wärmequellen in den Nutzraum kann es zu Kondensatbildung (z. B. an der Glastür) kommen.

Eingebrachte Energie	Beginn des Regelbereiches der Temperatur			
	allgemein	Beispiel: 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 + 13 °C	34 °C		
20 W	RT + 16 °C	37 °C		
*RT = Raumtemperatur				

Regalsystem



Abbildung 3-18. Komponenten des Regalsystems

Die Tragprofile (3/Abbildung 3-18) des Regalsystems sind in Abständen von 42 mm perforiert. Die Auflagebügel (2/Abbildung 3-18) können so für jede erforderliche Kulturgefäßgröße variabel eingesetzt werden. In die Einlagebleche (1/Abbildung 3-18) ist eine Kippsicherung und eine Auszugsbegrenzung eingearbeitet. Die Geräte **STERI-CYCLE i250** werden mit einem einteiligen unteren Einlageblech geliefert, siehe linke Seite der Abbildung 3-18, die Geräte **STERI-CYCLE i160** werden mit einem unteren Einlageblech mit Befüllausschnitt geliefert (4/Abbildung 3-18). Das Regalsystem wird im Kapitel "Inbetriebnahme" auf Seite 4-1 detailliert beschrieben.

Weitere Informationen zur Verwendung des Regalsystems mit 6-fach-Gasblende sind im Anhang zu finden.

Elektromechanisches Türverschluss-Kit

Das elektromechanische Türverschluss-Kit besteht aus einer Drehverriegelung (Detail A in Abbildung 3-19), die durch einen eingebauten Motor angetrieben wird, einem eingebauten mechanischen Notentriegelungshebel (Detail C) und einem Verschlusshaken (Detail B), der an die Türinnenseite montiert wird.

Das Türverschluss-Kit ist eine Sicherheitseinrichtung, die den Zugriff zum Nutzraum des Inkubators während der Steri-run-Dekontaminationsroutine verhindert. Diese Sicherheitseinrichtung verriegelt die Tür des Inkubators, sobald die Nutzraumtemperatur 65 °C erreicht/übersteigt. Bei Beendigung der Dekontaminationsroutine, entriegelt das Türverschluss-Kit die Türverriegelung, kurz nachdem die Temperatur unter 65 °C gesunken ist.



Abbildung 3-19. Türverriegelungshebel und Notentriegelung an der Unterseite des Inkubators

Bei Auftreten eines Stromausfalls bleibt der Hebel in seiner aktuellen Stellung. Nach Wiederherstellung der Stromversorgung und der Fortsetzung der Dekontaminationsroutine entriegelt der Hebel die Tür automatisch, sobald die Temperatur unter 65 °C sinkt. Die Tür kann ggf. jederzeit entriegelt werden, indem der Notentriegelungshebel (C) in Richtung Seitenwand gezogen wird. Der Notentriegelungshebel befindet sich an der Unterseite des Inkubators neben dem Netzschalter. Detail C in Abbildung 3-19 zeigt einen Ausschnitt durch die Seitenwand, um die Position des Hebels zu veranschaulichen.

Inbetriebnahme

Inhalt

- "Gerät akklimatisieren" auf Seite 4-2
- "Nutzraum aufbereiten" auf Seite 4-2
- "Füllstandsanzeiger "MAX" und Vorfilter einsetzen" auf Seite 4-3
- "Luftführung installieren" auf Seite 4-5
- "Regalsystem installieren" auf Seite 4-9
- "Geteilte Einlagebleche (optional) in STERI-CYCLE i250 einsetzen" auf Seite 4-11
- "Gas anschließen" auf Seite 4-11
- "Netzanschluss" auf Seite 4-16
- "USB-Schnittstelle anschließen:" auf Seite 4-17
- "Alarmkontakt anschließen:" auf Seite 4-18

Gerät akklimatisieren

VORSICHT Gerät akklimatisieren!



Vor Inbetriebnahme muss das Gerät akklimatisiert werden.

- Das Gerät vor dem Einschalten ca. 2 h im Betriebsraum bei voraussichtlicher Betriebsraumtemperatur aufstellen.
- Gerätetüren öffnen.

Nutzraum aufbereiten

 $\label{eq:CO2-Inkubator} Der\ CO_2 \mbox{-Inkubator} wird nicht in sterilem Zustand ausgeliefert. Vor der Aufnahme des Betriebes muss das Gerät dekontaminiert werden.$

Die folgenden Komponenten des Nutzraumes müssen dazu zuvor gereinigt und desinfiziert werden:

- Tragprofile
- Auflagebügel
- Vorfilter
- Luftführung (Airduct)
- Airbox
- Einlagebleche
- Oberflächen des Nutzraumes
- Glastürdichtung
- Glastür/Gasblende

Hinweis Dekontamination:

Detaillierte Maßnahmen zur Reinigung und Desinfektion des Gerätes sind in einem gesonderten Kapitel beschrieben (siehe "Reinigung und Desinfektion" auf Seite 8-1).

Füllstandsanzeiger "MAX" und Vorfilter einsetzen

Der Füllstandsanzeiger "MAX" und das Vorfilter lassen sich ohne Werkzeug einsetzen:

- 1. Sicherstellen, dass der Schlauch vom Befüll- und Ablassventil (3/Abbildung 4-2) abgezogen ist.
- 2. Prüfen, ob der Wasserablass (2/Abbildung 4-2) im vorderen Bereich des Wasserreservoirs frei ist; dieser wird für die Entleerung über das Befüll- und Ablassventil (3/Abbildung 4-2) an der Gerätevorderseite benötigt.
- Füllstandsanzeiger "MAX" (1/Abbildung 4-1) in den dafür vorgesehenen Schlitz in der Wasserreservoirabdeckung (5/Abbildung 4-1) einhängen.



Abbildung 4-1. Füllstandsanzeiger "MAX"



4. Vorfilter (4/Abbildung 4-2) in die Wasserreservoirabdeckung einsetzen.

Abbildung 4-2. Füllstandsanzeiger "MAX" und Vorfilter

Luftführung installieren

- 1. Oberteil (1/Abbildung 4-3) der Luftführung mit Rückteil (2/Abbildung 4-3) nach Abbildung 4-3, Schritte A-E zusammenfügen. Dabei darauf achten, dass die Positionierhilfslasche am Rückteil bei Schritt C in Abbildung 4-3 in das entsprechende Vierkantloch im Unterteil einrastet.
- 2. Steg am Fuß des Rückteils (2/Abbildung 4-3) auf die beiden Stehbolzen der Rückwand aufsetzen und Luftführung nach hinten kippen.
- 3. Die seitlichen Schlüssellöcher am Oberteil (Schritt G/Abbildung 4-3) in die Halteschrauben in der Nutzraumdecke einrasten.





Abbildung 4-3. Luftführung zusammenbauen



Abbildung 4-4 listet die Teile auf, aus denen die im Innenbehälter eingebaute Luftführung besteht.

Abbildung 4-4. Bauteile Luftführung STERI-CYCLE i160

Hinweis Bei den **Steri-CYCLE i160**-Ausführungen die Dichtungen der Luftführung nicht entfernen, bevor die Rückwand montiert ist. Die Luftführungen an der Rückwand sind wichtig für einen korrekten Luftstrom in den **Steri-CYCLE i160**-Ausführungen mit einem Innenbehälter aus Edelstahl.

HEPA-Filter und Wasserreservoirabdeckung einbauen

Das HEPA-Filter (2/Abbildung 4-5) wird von unten in die Airbox (1/Abbildung 4-5) eingesetzt. Die Airbox sitzt auf einem Sockel auf der Wasserreservoirabdeckung (2/Abbildung 4-6) und wird an den Lüftereinlass geschoben.



Abbildung 4-5. HEPA-Filter und Airbox zusammenbauen

- 1. HEPA-Filter (2/Abbildung 4-5) auf gerade Oberfläche legen.
- 2. Airbox (1/Abbildung 4-5) nach links ankippen und mit den Laschen auf der linken Seite (3/Abbildung 4-5) in die entsprechenden Nuten im HEPA-Filter (4/Abbildung 4-5) einführen.
- 3. Laschen (5/Abbildung 4-5) auf der rechten Seite der Airbox in die Rasten (6/Abbildung 4-5) des HEPA-Filters einrasten lassen.
- 4. Dichtung (8/Abbildung 4-5) in Nut (7/Abbildung 4-5) an Rohrstutzen der Airbox einsetzen und rundherum festdrücken.

5. Airbox (1/Abbildung 4-6) auf Sockel (2/Abbildung 4-6) der Wasserreservoirabdeckung aufsetzen.



Abbildung 4-6. Airbox auf die Wasserreservoirabdeckung aufsetzen

- 6. Wasserreservoirabdeckung auf Nutzraumboden aufsetzen.
- 7. Die Wasserreservoirabdeckung an der Frontseite anheben und in Richtung Rückwand schieben (1/Abbildung 4-7).



Abbildung 4-7. Airbox einbauen

- 8. Wasserreservoirabdeckung bis zum Anschlag in Richtung Rückwand schieben. Die Abdeckung gleitet in ihre Endposition im Reservoir und der Rohrstutzen der Airbox in den Lüfterauslass.
- 9. Vorderkante der Wasserreservoirabdeckung in das Wasserreservoir gleiten lassen (2/Abbildung 4-7). Dabei gleitet der Stutzen der Airbox in den Lüfterauslass.

Regalsystem installieren



Abbildung 4-8. Ein-/Ausbau Regalsystem

Zur Installation des Regalsystems wird kein Werkzeug benötigt. Die Tragprofile werden durch Federdruck gehalten. Die Auflagebügel werden in die Tragprofile eingehängt, die Einlagebleche auf die Auflagebügel geschoben.

Tragprofile ein-/ausbauen

Die Tragprofile werden durch die Prägungen seitlich geführt und fixiert. Die Haltefedern der Tragprofile müssen dabei nach oben zeigen.

- 1. Tragprofil auf die untere Prägung setzen und an die Seitenwand des Nutzraumes klappen, so dass das Tragprofil über den beiden Prägungen sitzt.
- 2. Die Haltefeder hinter die obere Prägung klemmen.
- 3. Zum Ausbau der Tragprofile die Haltefeder an der Lasche nach unten aus der Prägung ziehen und das Tragprofil herausnehmen.

Auflagebügel einsetzen



Abbildung 4-9. Auflagebügel einsetzen

- 1. Auflagebügel in die Perforation des Tragprofils stecken, so dass der Auflagestab nach unten zeigt.
- 2. Sicherstellen, dass die beiden Vertikalstücke des Auflagebügels am Tragprofil anliegen.

Einlagebleche einschieben:

- 1. Die Einlagebleche mit der Kippsicherung (5/Abbildung 4-9) zur Geräterückwand zeigend auf den Auflagebügel schieben. Die Kippsicherung dient gleichzeitig als Führung des Einlagebleches.
- 2. Einlageblech leicht anheben, so dass die Auszugsbegrenzung (6/Abbildung 4-9) über den Auflagebügel geführt werden kann.
- 3. Sicherstellen, dass der Auflagebügel frei in den beiden Kippsicherungen läuft.

Hinweis Beim **STERI-CYCLE i160**-Inkubator das Einlageblech mit Befüllvorrichtung an unterster Stelle einschieben. Bei der 6-fach-Gasblende werden die Auflagebügel nur in die Vierkantlöcher eingesetzt, siehe Anhang.

Gerät nivellieren

- 1. Eine Wasserwaage auf das mittlere Einlageblech bzw. auf die Rollenaufnahme legen.
- Die verstellbaren Gerätefüße mit dem mitgelieferten Schraubenschlüssel (Schlüsselweite 24 mm) verdrehen, so dass das Einlageblech nach allen Richtungen waagerecht ausgerichtet ist. Die Höheneinstellung der Gerätefüße sollte von links nach rechts und von hinten nach vorne erfolgen.

Geteilte Einlagebleche (optional) in STERI-CYCLE i250 einsetzen



Abbildung 4-10. Einbau geteilter Einlagebleche

Ist **STERI-CYCLE i250** mit der optionalen 6-fach-Gasblende und den 6-fach geteilten Einlageblechen ausgestattet, werden statt der einteiligen Horden der Normalausstattung drei Tragrahmen mit jeweils zwei Einlageblechen auf die Auflagebügel der seitliche Tragprofile aufgesetzt.

- 1. Beide Einlagebleche (1/Abbildung 4-10) auf den Tragrahmen (2/Abbildung 4-10) aufsetzen (Schritt A).
- 2. Tragrahmen (2/Abbildung 4-10) in die Auflagebügel (3/Abbildung 4-10) einhängen (Schritt B).

Gas anschließen

Hinweis Qualität des Gases:

Die Gase müssen eines der folgenden Qualitätsmerkmale aufweisen:

- Reinheit min. 99,5 %
- medizinische Gasqualität.

VORSICHT Überdruck!



Gas darf am Gerät mit maximal 1 bar Betriebsdruck angelegt werden. Wird das Gas mit einem höheren Druck eingeleitet, besteht die Gefahr, dass die geräteinternen Ventile nicht richtig schließen und die Regelung der Gaszufuhr deshalb nicht einwandfrei funktioniert. Die Gaszufuhr auf einen Bereich von min. 0,8 bar und max. 1 bar einstellen und sicherstellen, dass dieser Vordruck nicht verändert werden kann!

Gasdruckschläuche montieren



Abbildung 4-11. Gasdruckschläuche montieren

Die Gaszuleitung von der Gasversorgungsanlage zum Gerät wird durch die mitgelieferten flexiblen Gasdruckschläuche hergestellt:

- 1. Gasdruckschlauch auf den Anschlussstutzen der Gasversorgungsanlage stecken.
- 2. Schutzkappe (3/Abbildung 4-11) des Gaseinlassfilters abnehmen.
- Schlauchschelle (1/Abbildung 4-11) auf den Gasdruckschlauch (2/Abbildung 4-11) schieben und den Gasdruckschlauch auf den Anschlussstutzen (4/Abbildung 4-11) des Gaseinlassfilters (5/Abbildung 4-11) stecken.
- 4. Gasdruckschlauch mit der Schlauchschelle am Anschlussstutzen des Gaseinlassfilters fixieren.

VORSICHT Druckausgleichsöffnung

Um einen ständigen Druckausgleich zu ermöglichen, darf die Druckausgleichsöffnung nicht an ein Abluftsystem angeschlossen sein. Das Rohr der Druckausgleichsöffnung darf weder verlängert noch umgeleitet werden.


Gasanschluss ohne Gasmonitoring

Abbildung 4-12. Gasanschluss ohne Gasmonitoring

Die Gaszuleitung von der Gasversorgungsanlage zum Gerät wird durch die mitgelieferten flexiblen Gasdruckschläuche hergestellt:

CO₂-Anschluss:

• Bei einem Gerät mit CO₂-Anschluss wird die Gasversorgung an das Gaseinlassfilter (2/Abbildung 4-12) gemäß Anschlussschema (1/Abbildung 4-12) angeschlossen.

Kombinierter CO₂- und O₂-/N₂-Anschluss (optional):

Bei einem kombinierten CO_2 - $/O_2$ - $/N_2$ -Anschluss werden die Zuleitungen gemäß Anschlussschema (3/Abbildung 4-12) folgendermaßen verlegt:

• O₂-/N₂-Versorgung an das Gaseinlassfilter oben (4/Abbildung 4-12),

• die CO₂-Versorgung an das Gaseinlassfilter unten (5/Abbildung 4-12).

Hinweis Manuelle Gasüberwachung:

Bei Geräten ohne Gasmonitoring (optional) findet keine automatische Gasüberwachung statt. Der Füllstand der Gasflaschen muss deshalb täglich überprüft werden.

CO₂-Anschluss mit Gasmonitoring-System (optional)



Abbildung 4-13. CO₂-Anschluss mit Gasmonitoring-System (optional)

Die Gaszuleitung von der Gasversorgungsanlage zum Gerät wird durch die mitgelieferten flexiblen Gasdruckschläuche hergestellt.

Geräte mit CO_2 -Anschluss, die mit dem optionalen Gasmonitoring-System ausgerüstet sind, werden gemäß Anschlussschema (7/Abbildung 4-13) angeschlossen.

CO₂-Anschluss:

Bei einem Gerät, das mit dem Prozessgas CO₂ betrieben wird und das mit dem optionalen Gasmonitoring-System ausgerüstet ist, wird die Gasversorgung folgendermaßen angeschlossen:

- Position 1/Abbildung 4-13: Verteileranschluss f
 ür die CO₂-Gasversorgung von bis zu drei weiteren Ger
 äten. Zur Weiterschleifung der Gasversorgung die Ger
 äte durch einen Gasdruckschlauch miteinander verbinden.
- Position 2/Abbildung 4-13: Die Gasversorgung von Druckgasbehälter B am oberen Anschluss des Gasmonitoring-Systems anlegen.
- Position 3/Abbildung 4-13: Die Gasversorgung von Druckgasbehälter A am unteren Anschluss des Gasmonitoring-Systems anlegen.

 Position 4 & 5/Abbildung 4-13: Der Ausgang des Gasmonitoring-Systems ist werkseitig bereits durch einen kurzen Gasdruckschlauch (Position 6/Abbildung 4-13) mit dem Gaseinlassfilter verbunden.

Hinweis Weiterschleifung der Gasversorgung: Durch Weiterschleifung der Gaszufuhr können insgesamt maximal 4 Geräte durch je eine CO₂-Gasversorgung versorgt werden.

Kombinierter CO_2 - und O_2/N_2 -Anschluss mit Gasmonitoring-System (optional)



Abbildung 4-14. Kombinierter CO₂- und O₂-/N₂-Anschluss mit Gasmonitoring-System (optional)

Die Gaszuleitung von der Gasversorgungsanlage zum Gerät wird durch die mitgelieferten flexiblen Gasdruckschläuche hergestellt.

Geräte, die mit dem kombinierten CO_2 - $/O_2$ - $/N_2$ -Anschluss und mit dem optionalen Gasmonitoring-System ausgerüstet sind, werden gemäß Anschlussschema (10/Abbildung 4-14) angeschlossen.

O₂-/N₂-Versorgung:

- Verteileranschluss f
 ür die O₂-/N₂-Gasversorgung von bis zu drei weiteren Ger
 äten (verdeckt in Abbildung 4-14). Zur Weiterschleifung der Gasversorgung die Ger
 äte durch einen Gasdruckschlauch miteinander verbinden.
- Position 2/Abbildung 4-14: Die Gasversorgung von Druckgasbehälter B am Anschluss des Gasmonitoring-Systems anlegen.
- Position 3/Abbildung 4-14: Die Gasversorgung von Druckgasbehälter A am Anschluss des Gasmonitoring-Systems anzulegen.
- Position 1/Abbildung 4-14: Der Ausgang des Gasmonitoring-Systems ist werkseitig bereits durch einen kurzen Gasdruckschlauch mit dem Gaseinlassfilter (Position 8/Abbildung 4-14) verbunden.

CO₂-Versorgung:

- Position 4/Abbildung 4-14: Verteileranschluss für die CO₂-Gasversorgung von bis zu drei weiteren Geräten. Zur Weiterschleifung der Gasversorgung die Geräte durch einen Gasdruckschlauch miteinander verbinden.
- Position 5/Abbildung 4-14: Die Gasversorgung von Druckgasbehälter B am Anschluss des Gasmonitoring-Systems anlegen.
- Position 6/Abbildung 4-14: Die Gasversorgung von Druckgasbehälter A am Anschluss des Gasmonitoring-Systems anlegen.
- Position 7/Abbildung 4-14: Der Ausgang des Gasmonitoring-Systems ist werkseitig bereits durch einen kurzen Gasdruckschlauch mit dem Gaseinlassfilter verbunden.

Hinweis Weiterschleifung der Gasversorgung:

Durch Weiterschleifung der Gaszufuhr können insgesamt maximal 4 Geräte (unabhängig vom Gerätetyp) durch je eine CO₂-Gasversorgung bzw. O₂-/N₂-Gasversorgung versorgt werden.

Netzanschluss



Abbildung 4-15. Netzanschluss



WARNUNG Stromschlag!

Die Berührung stromführender Teile kann zu einem lebensgefährlichen Stromschlag führen. Stecker und Stromkabel vor dem Netzanschluss auf Beschädigung überprüfen. Beschädigte Anschlusskomponenten dürfen nicht zum Netzanschluss verwendet werden!

Das Gerät an ein vorschriftsmäßig installiertes und geerdetes Stromnetz mit Absicherung durch einen Leitungsschutzschalter B 16 anschließen.

Netzanschluss herstellen:

- Überprüfen Sie vor dem Anschluss an das Netz, ob die Spannungswerte der Steckdose mit den Angaben auf dem Typenschild an der Seite links vom Netzschalter des Gerätes übereinstimmen. Stimmen die Angaben für Spannung (V) und maximalen Strom (A) nicht überein, darf das Gerät nicht angeschlossen werden.
- 2. Den Kaltgerätestecker in die Steckdose am Schaltkasten des Gerätes stecken.
- 3. Den Schutzkontaktstecker des Netzkabels in eine ordnungsgemäß geerdete und abgesicherte Steckdose stecken.
- 4. Stellen Sie sicher, dass keine Zug- oder Druckkräfte auf das Netzkabel einwirken.
- 5. Installation des Netzanschlusses: Der Netzstecker muss für den Benutzer klar erkennbar und jederzeit frei zugänglich sein. Der Stecker der Netzzuleitung bildet die allpolige Trennvorrichtung.

USB-Schnittstelle anschließen:

Die Geräte sind serienmäßig mit einer USB-Schnittstelle ausgerüstet. Die Verbindung zum PC wird über ein handelsübliches USB 1- oder USB 2-Kabel (Inkubatorseite: USB-Stecker Typ B, PC-Seite: USB-Stecker Typ A) hergestellt.

Die USB-Schnittstelle entspricht dem Standard USB 1.1 und ist mit den Standards USB 2.0 und 3.0 (full speed) kompatibel. Zur Datenkommunikation zwischen PC und Inkubator über die USB-Schnittstelle muss auf dem PC ein entsprechender Treiber installiert werden, sofern dieser nicht von der aktuell installierten Windows-Version bereitgestellt wird. Die Installation des Treibers ist im "Datenkommunikation" auf Seite 12-1 beschrieben.

Die Übertragungsgeschwindigkeit der Schnittstelle kann bei Bedarf innerhalb der definierten Baudraten (9.600, 19.200, 38.400, 57.600 Baud) verändert werden. Die Einstellung der Baudrate ist im Abschnitt "Baudrate USB-Schnittstelle einstellen" auf Seite 6-23 beschrieben.

Alarmkontakt anschließen:

Hinweis Facharbeiten:

Thermo Scientific gewährleistet die Sicherheit und Funktionstüchtigkeit des Gerätes nur, wenn Installationen und Instandsetzungsarbeiten fachgerecht ausgeführt werden. Der Anschluss des Gerätes an ein externes Alarmsystem darf nur von ausgebildetem und autorisiertem Fachpersonal der Elektrotechnik/Fernmeldetechnik ausgeführt werden!

Funktion:

Beim Auftreten von Systemfehlern und Fehlern in den Regelkreisen von Temperatur oder Gas wird an das angeschlossene Melde-/Überwachungssystem eine Alarmmeldung abgegeben. Der potentialfreie Kontakt (1 Wechsler) ist für folgende Stromkreise dimensioniert:

Alarmrelais:

Stromkreis	Spannung	Externe Absicherung
Stromkreise mit Netzspannung	max. 250 V ~	max. 6 A
SELV – Stromkreise	25 V ~	max. 2 A
(vgl. VDE 0100, Teil 410)	60 V =	max. 1 A
SELV E – Stromkreise	50 V ~	max. 1 A
(vgl. VDE 0100, Teil 410)	120 V =	max. 0,5 A
Detrick envetive de	Kontokt 2 0	Kenteld 0 4
Betriebszustande	Kontakt 3 - 2	Kontakt 3 - 1
Betriebszustande Betriebszustand Netzüberwachung "aus"	X	C
Betriebszustande Netzüberwachung "aus" Betriebszustand	коптакт 3 - 2 Х О	коптакт 3 - 1 О Х
Betriebszustande Netzüberwachung "aus" Betriebszustand Netzüberwachung "ein"	коптакт 3 - 2 Х О	CONTART 3 - 1 O X
Betriebszustande Betriebszustand Netzüberwachung "aus" Betriebszustand Netzüberwachung "ein" Fehler	X 0 O 0	Kontakt 3 - 1 O X X
Betriebszustande Betriebszustand Netzüberwachung "aus" Betriebszustand Netzüberwachung "ein" Fehler Netzüberwachung "aus"	X 0 O 0	Kontakt 3 - 1 O X X
Betriebszustande Betriebszustand Netzüberwachung "aus" Betriebszustand Netzüberwachung "ein" Fehler Netzüberwachung "aus" Fehler	X 0 O X X X	Kontakt 3 - 1 O X X O
Betriebszustande Betriebszustand Netzüberwachung "aus" Betriebszustand Netzüberwachung "ein" Fehler Netzüberwachung "aus"	X O O X	Kontakt 3 - 1 O X X O

Hinweis Schaltstruktur:

Das Alarmrelais schaltet bei allen, von den Regelkreisen gemeldeten Fehlern (siehe "Fehlermeldungen" auf Seite 6-45).

Anschlussbeispiel:



Abbildung 4-16. Beispiel Anschluss Alarmkontakt

Der Stecker zum Anschluss des Verbindungskabels gehört nicht zum Lieferumfang, kann jedoch separat bestellt werden. Die Werte für die Betriebsspannung und Absicherung der externen Stromkreise des Meldesystems sind in der Tabelle beschrieben.

- 1. Die einzelnen Litzen des Verbindungskabels gemäß den Zuordnungen im Stromanschlussplan anklemmen.
- 2. Stecker des Verbindungskabels zum externen Meldesystem in die Schnittstelle am Schaltkasten an der Geräterückseite stecken.

Schnittstelle 4-20mA anschließen:

Die folgende Abbildung zeigt die Belegung der Messdatenschnittstelle 4-20mA:

Abbildung 4-17. Anschlussbelegung der Schnittstelle 4-20 mA

Hinweis Facharbeiten:

Thermo Scientific gewährleistet die Sicherheit und Funktionstüchtigkeit des Gerätes nur, wenn Installationen und Instandsetzungsarbeiten fachgerecht ausgeführt werden.

Der Anschluss des Gerätes an ein externes Alarmsystem darf nur von ausgebildetem und autorisiertem Fachpersonal der Elektrotechnik/Fernmeldetechnik ausgeführt werden!

Die Nachrüstung der Messdatenschnittstelle 4-20 mA darf ausschließlich durch den Technischen Service von Thermo Fisher Scientific vorgenommen werden.

Betrieb

Inhalt

- "Gerät vorbereiten" auf Seite 5-1
- "Betrieb aufnehmen" auf Seite 5-2

Gerät vorbereiten

Das Gerät darf nur für den laufenden Betrieb freigegeben werden, wenn alle wichtigen Maßnahmen zur Inbetriebnahme (siehe "Inbetriebnahme" auf Seite 4-1) durchgeführt wurden.

Geräte-Check:

Bevor der Betrieb aufgenommen wird, muss in einem Geräte-Check der Zustand folgender Bauteile geprüft werden:

- Gasschläuche müssen dicht auf den Anschlussfiltern sitzen und mit Schlauchschellen gesichert sein.
- Die Rohrdurchführung muss verschlossen sein.
- Druckausgleichsöffnung muss durchlässig sein, der Einsatz der Druckausgleichsöffnung im Nutzraum installiert sein.
- Die Dichtung der Glastür darf nicht beschädigt sein.
- Die Messöffnung der Glastür/Gasblende muss verschlossen sein.
- Die Komponenten des Regalsystems müssen sicher eingebaut sein.
- Die Airbox mit HEPA-Filter und die Luftführung müssen ordnungsgemäß eingebaut sein.

Nutzraum des Gerätes dekontaminieren:



VORSICHT Betriebstemperatur-Grenzwerte des Cell Lockers! Vor der Durchführung eines Dekontaminationslaufs sind die Cell Locker zu entfernen. Die max. Betriebstemperatur beträgt 121°C/250 °F.

• Steri-run-Dekontaminationsroutine durchführen (siehe "Steri-run-Dekontaminationsroutine" auf Seite 8-7) oder den Nutzraum entsprechend den vom Betreiber festgelegten Hygienerichtlinien dekontaminieren.

Hinweis Hygienerichtlinien: Zum Schutz der Kulturen muss der Nutzraum des Gerätes entsprechend den vom Betreiber festzulegenden Hygienerichtlinien vor jedem Betriebseinsatz gereinigt und desinfiziert werden.

Wasservorrat: siehe "Relative Feuchte:" auf Seite 3-6.

Bei Unterschreiten der Mindestfüllmenge muss während des Arbeitsprozesses der Wasservorrat ergänzt werden können.

Füllmenge STERI-CYCLE i160 und STERI-CYCLE i250: 31

Betrieb aufnehmen

- 1. Glastür oder Gasblende öffnen.
- 2. Vorfilter (2/Abbildung 5-1) herausnehmen.
- Einlageblech mit Befüllausschnitt (STERI-CYCLE i160): Klappe öffnen (3/Abbildung 5-1). Einteiliges Einlageblech (STERI-CYCLE i250): Unteres Einlageblech herausnehmen, wenn der Platz für das Gefäß zum Befüllen nicht ausreicht.
- 4. Ausreichend aufbereitetes Wasser in die Bodenwanne (unter Abdeckung Position 1 in Abbildung 5-1) des Nutzraumes gießen.



Abbildung 5-1. Wasserreservoir

Vorgehensweise beim Befüllen mit Wasser



Abbildung 5-2. Befüll- und Ablassventil des Inkubators

Der Inkubator kann mit einem Befüllzylinder über das Wasserventil erneut mit Wasser befüllt werden.



Abbildung 5-3. Befüllzylinder

Die Haken der Befüllvorrichtung in die Verschlüsse der Glastüren einhängen (siehe oben). Den Schlauch an das Befüll- und Ablassventil anschließen.

Der Befüllzylinder bestimmt nach dem Prinzip der kommunizierenden Röhren die Höhe des Wasserstands im Inkubator.

Die minimale Füllmenge beträgt 0,5 l, die maximale Füllmenge beträgt 3,5 l. Der Zylinder enthält ca. 0,4 l Wasser.

- Den Zylinder in den Wasserablass einsetzen. Einige Zeit warten, bis sich der Wasserstand stabilisiert hat, sodass dieser abgelesen werden kann.
- Um das Befüllen durch die Einwirkung der Schwerkraft zu ermöglichen, Zylinder bis zum Maximum mit sterilem destilliertem Wasser befüllen. Dieser Vorgang muss ggf. mehrmals wiederholt werden.
- Wasser hinzufügen, bis die maximale Füllstandsanzeige erreicht ist.

5. Die Füllmenge darf die maximale Füllstandsmarke "MAX" (1/Abbildung 5-1) nicht überschreiten. Als maximale Füllmenge für das Wasserreservoir (5/Abbildung 5-1) sind 3 l zulässig.



Abbildung 5-1. Füllstandsanzeiger "MAX"

- 6. Überschüssiges Wasser vom Deckel des Wasserreservoirs abwischen.
- 7. Vorfilter (2/Abbildung 5-1) wieder einsetzen.
- 8. Unteres Einlageblech wieder einsetzen bzw. Klappe des Befüllausschnitts schließen.
- 9. Sicherstellen, dass die Ventile der CO₂-/O₂-/N₂-Versorgungsanlage geöffnet sind.
- 10. Gerät am Netzschalter einschalten.
- 11. Sollwerte für Temperatur und CO₂-/O₂-Gehalt am iCan[™] Touchscreen einstellen.

Gerät starten:

- 12. Gerät mit auto-start starten (siehe "auto-start aktivieren" auf Seite 6-14).
- 13. Die Verlaufsanzeige auto-start wird am Display eingeblendet, die automatische Startroutine läuft.
- 14. Die Temperaturregelung regelt auf den eingestellten Temperatur-Sollwert, die Feuchte wird aufgebaut.
- 15. Nachdem Temperatur und relative Feuchte konstant sind, erfolgt der automatische Abgleich des CO₂-/O₂-Messsystems.
- 16. Die CO₂-/O₂-Regelung begast auf den eingestellten CO₂-/O₂-Sollwert.
- 17. Ist die Routine auto-start beendet, wird die Verlaufsanzeige ausgeblendet und das Hauptmenü angezeigt. Das Gerät ist betriebsbereit.

Gerät beschicken:

18. Nutzraum mit Kulturen beschicken.

Hinweis

Dauer der auto-start-Routine:

Bei kaltem Gerät und niedrigen Umgebungstemperaturen kann die auto-start-Routine bis zu 10 Stunden dauern.

Beschickung:

Damit eine ausreichende Luftzirkulation und eine gleichmäßige Erwärmung der Proben möglich ist, sollte die Beschickungsfläche im Nutzraum max. zu 70 % genutzt werden. Großflächige Gegenstände oder Geräte mit Wärmeabgabe im Nutzraum können die Wärmeverteilung beeinträchtigen. Großflächige Gegenstände oder Geräte mit Wärmeabgabe im Nutzraum können die Wärmeverteilung beeinträchtigen.

5 Betrieb Vorgehensweise beim Befüllen mit Wasser

Bedienung

Inhalt

- "Netzschalter" auf Seite 6-2
- "Bedienfeld und Bedienstruktur" auf Seite 6-3
- "Werkseinstellungen der Regler des iCan™ Touchscreens" auf Seite 6-7
- "Aufwärmphase der Sensoren der Regelkreise" auf Seite 6-7
- "Verhalten der Tasten bei Einstellungen" auf Seite 6-8
- "Temperatur-Sollwert einstellen" auf Seite 6-8
- "CO2-Sollwert einstellen" auf Seite 6-9
- "O2-Sollwert einstellen" auf Seite 6-10
- "auto-start-Funktion" auf Seite 6-12
- "Steri-run aufrufen" auf Seite 6-16
- "Benutzerkonfiguration" auf Seite 6-17
- "Verlaufsanzeige skalieren" auf Seite 6-43
- "Fehlermeldungen" auf Seite 6-45
- "Maßnahmen nach einem Stromausfall" auf Seite 6-47

Netzschalter



Abbildung 6-1. Netzschalter

Der Netzschalter ist in die Geräteseitenwand eingelassen.

- Gerät einschalten: Netzschalter drücken, der Schalter wird beleuchtet.
 - Nach einem kurzen Signalton und einer kurzen Dunkelphase wird das Display des Tastbildschirms angezeigt.
 - Die Sensoren der Regelkreise durchlaufen die Aufwärmphase ("Aufwärmphase der Sensoren der Regelkreise" auf Seite 6-7).
- Gerät ausschalten: Netzschalter drücken, die Schalteranzeige erlischt.

Bedienfeld und Bedienstruktur

Das Bedienfeld ist ein Tastbildschirm (iCan™ Touchscreen) und kann durch leichten Druck mit einem Finger oder einem stumpfen Stift bedient werden.



Abbildung 6-2. Hauptbildschirm: Tastsensitive Bildschirmbereiche

Folgende Bildschirmbereiche des Bedienfelds sind druckempfindlich und können bedient werden:

- Temperatur-Anzeigefeld T,
- CO₂-Anzeigefeld CO₂,
- Anzeigefeld Wasserstand RH,
- O₂-Anzeigefeld (optional),
- Icon-Leiste mit Symbolen zur Anzeige von Betriebszuständen und zum Direktzugriff auf installierte Optionen (siehe auch "Iconerklärung" auf Seite 6-40),
- steri-run-Taste,
- Menü-Taste,
- auto-start-Taste.

Hinweis Erweiterter druckempfindlicher Bereich: Zur Bestätigung einer Fehlermeldung kann der gesamte Tastbildschirm als druckempfindlicher Bereich benutzt werden.

Ausstattung ohne O₂-/N₂-Regelung:



Abbildung 6-3. iCan[™] Touchscreen ohne Gasversorgung O₂/N₂

Funktionstasten und Istwertanzeigen des Bedienfeldes bei einer Geräte
ausstattung ohne Gasversorgung $\rm O_2/N_2\!\!:$

- 1. Anzeigefeld Arbeitsraumtemperatur mit Istwert (große Zahl in der Mitte), Sollwert (kleine Zahl rechts oben) und physikalischer Einheit (rechts unten)
- 2. Anzeigefeld CO²-Konzentration mit Istwert (große Zahl in der Mitte), Sollwert (kleine Zahl rechts oben) und physikalischer Einheit (rechts unten)
- 3. Aktivitätsanzeige HEPA-Filter (Symbole für weitere Optionen siehe "Iconerklärung" auf Seite 6-40)
- 4. Taste zum Start der Steri-run-Dekontaminationsroutine
- 5. Taste zum Öffnen der Menü-Navigation
- 6. Taste zum Start der auto-start-Routine
- 7. Anzeigefeld Wasserstand RH



Ausstattung mit kombinierter CO₂-/O₂-/N₂-Regelung (optional):

Abbildung 6-4. iCan™ Touchscreen mit kombiniertem Gasanschluss

Funktionstasten und Istwertanzeigen des Bedienfeldes bei einer Geräteausstattung mit kombinierter Gasversorgung CO₂/O₂/N₂:

- 1. Anzeigefeld Arbeitsraumtemperatur mit Istwert (große Zahl in der Mitte), Sollwert (kleine Zahl rechts oben) und Einheit (rechts unten)
- 2. Anzeigefeld CO₂-Konzentration mit Istwert (große Zahl in der Mitte), Sollwert (kleine Zahl rechts oben) und Einheit (rechts unten)
- 3. Aktivitätsanzeige HEPA-Filter (Symbole für weitere Optionen siehe "Iconerklärung" auf Seite 6-40)
- 4. Taste zum Start der Steri-run-Dekontaminationsroutine
- 5. Taste zum Öffnen der Menü-Navigation
- 6. Taste zum Start der auto-start-Routine
- 7. Anzeigefeld O₂-Konzentration mit Istwert (große Zahl in der Mitte), Sollwert (kleine Zahl rechts oben) und physikalischer Einheit (rechts unten)
- 8. Anzeigefeld Wasserstand RH

Gliederung der Bedienebenen

Die Bedienung ist in drei Ebenen gegliedert:

- A: Direktzugriff zu den Einstellungen der Regelkreise: Temperatur-, CO₂-, O₂-Sollwert,
- B: Start der Geräteroutinen steri-run bzw. auto-start,
- C: Navigation durch die Untermenüs zur Konfiguration des Gerätes.



Abbildung 6-5. Übersicht Menüstruktur

Werkseinstellungen der Regler des iCan™ Touchscreens

Die Geräte sind bei Auslieferung auf folgende Sollwerte eingestellt:

- Temperatur: 37 °C
- CO₂-Gehalt: 5,0 %
- O₂-Gehalt (optional): 21,0 %

Hinweis CO₂-/O₂-Regelung:

Da die CO₂-Konzentration der Luft nahezu 0 % beträgt, sind die CO₂-Regelung sowie die Fehlerüberwachung des Regelkreises bei einem Sollwert von 0 % deaktiviert. Da die O₂-Konzentration der Luft 21 % beträgt, ist die O₂-Regelung, sowie die Fehlerüberwachung des Regelkreises, bei einem Sollwert von 21 % deaktiviert.

Aufwärmphase der Sensoren der Regelkreise

Nach dem Einschalten des Gerätes durchlaufen die Sensoren der Regelkreise während des Startvorgangs eine Aufwärmphase mit unterschiedlicher Dauer:

- 1. Regelkreis Temperatur: 10 s
- 2. Aufwärmzeit CO2-Sensor (WLD und IR): 5 Min.
- 3. O₂-Regelkreis: 5 Min.



Abbildung 6-6. Anzeige Aufwärmphase

Der Startvorgang wird durch einen Signalton angekündigt. Während der Aufwärmphase werden Punkte (...) anstelle von Werten in den Anzeigefeldern angezeigt:

- Temperatur-Anzeigefeld,
- CO₂-Anzeigefeld und
- O₂-Anzeigefeld

Nach durchlaufener Aufwärmphase werden die Istwerte der Regelkreise angezeigt.

Hinweis CO₂-Begasung:

Während der 5-minütigen Aufwärmphase des O₂-Regelkreises erfolgt keine CO_2 -Begasung des Nutzraumes und keine Überwachung des $_{CO}$ 2-Regelkreises.

Verhalten der Tasten bei Einstellungen

Ein Wert kann durch Tastendruck stufenweise erhöht oder vermindert werden:

- Durch Dauerdruck auf die jeweilige Taste oder Taste + kann in einen Schnelldurchlauf umgeschaltet werden,
- nach ca. 3 s Dauerdruck wird die Geschwindigkeit des Schnelldurchlaufes zusätzlich erhöht.

Hinweis Speichern der Einstellungen: Geänderte Werte werden gespeichert, wenn sie durch Drücken der Taste **Enter** bestätigt werden.

Rücksetzung der Einstellungen:

Erfolgt 30 s lang keine Benutzeraktion (Berührung der sensitiven Bereiche und Tasten), wird das Menü automatisch verlassen und die zuletzt bestätigte Einstellung wird übernommen.

Temperatur-Sollwert einstellen

- 1. Taste **Temperatur** drücken.
- Das Temperatur-Menü (Abbildung 6-7) wird angezeigt.





Abbildung 6-7. Temperatur-Anzeigefeld und Temperatur-Auswahlmenü

- 2. Temperatur-Menü wieder verlassen:
- Taste Ende drücken.
- 3. Temperatur-Sollwert einstellen:
- Taste **Sollwert** drücken.





Sollwert erhöhen:

• Taste + drücken.

Sollwert verringern:

- Taste drücken.
- 4. Sollwert übernehmen und speichern:
- Taste Enter drücken.
- Es erfolgt ein Rücksprung ins Hauptmenü. Der aktuelle im Nutzraum gemessene Istwert wird am Temperatur-Anzeigefeld angezeigt.

CO₂-Sollwert einstellen

- 1. Taste CO₂-Anzeigefeld drücken.
- Das CO₂-Menü wird angezeigt.



Abbildung 6-9. CO₂-Anzeigefeld und CO₂-Menü

- 2. CO₂-Menü wieder verlassen:
- Taste Ende drücken.

- 3. CO₂-Sollwert einstellen:
- Taste **Sollwert** drücken.

Sollwert erhöhen:

• Taste + drücken.

Sollwert verringern:

- Taste drücken.
- 4. Sollwert übernehmen und speichern:
- Taste Enter drücken.
- Es erfolgt ein Rücksprung ins Hauptmenü. Der aktuelle, im Nutzraum gemessene Istwert wird am CO₂-Anzeigefeld angezeigt.



Abbildung 6-10. CO₂-Sollwert einstellen

Hinweis CO₂-Regelkreis deaktivieren:

Zur Deaktivierung der CO₂-Regelung wird der Soll-Wert auf 0 % gesetzt. Ist der Regelkreis deaktiviert, wird auch keine Fehlerüberwachung durchgeführt.

Gasmonitoring:

Der Schaltzustand des CO₂-Regelkreises hat keinen Einfluss auf die Funktion des optionalen Gasmonitorings. Auch bei ausgeschalteter CO₂-Regelung ist das Gasmonitor-System aktiv.

Hinweis

Den Innenbehälter nach der Änderung des CO₂-Sollwertes belüften, damit keine Alarmmeldung auftritt.

O₂-Sollwert einstellen

Diese Einstellung ist nur bei Ausstattung mit optionaler O₂-/N₂-Regelung möglich.

- 1. Taste O2-Anzeigefeld drücken.
- Das O₂-Menü wird angezeigt.



Abbildung 6-11. O2-Anzeigefeld und O2-Menü

- 2. O₂-Menü verlassen:
- Taste Ende drücken.
- 3. O₂-Sollwert einstellen:
- Taste **Sollwert** drücken.

Sollwert erhöhen:

• Taste + drücken.

Sollwert verringern:

- Taste drücken.
- 4. Sollwert übernehmen und speichern:
- Taste **Enter** drücken.
- Es erfolgt ein Rücksprung ins Hauptmenü. Der aktuelle, im Nutzraum gemessene Istwert wird am O₂-Anzeigefeld angezeigt.



Abbildung 6-12. O₂-Sollwert einstellen

Hinweis

Den Innenbehälter nach der Änderung des O₂-Sollwertes belüften, damit keine Alarmmeldung auftritt.

Hinweis Werkseinstellungen:

Abhängig vom Typ des O₂-Sensors ist werkseitig einer der beiden O₂-Regelbereiche voreingestellt: Regelbereich I: 1 % - 21 % Regelbereich II: 5 % - 90 %

Verwendung der Prozessgase:

Bei Sollwerten unter 21 % O_2 ist das Gerät an eine Stickstoffversorgung anzuschließen. Bei Sollwerten über 21 % O_2 ist das Gerät an eine Sauerstoffversorgung anzuschließen. Bei einem Sollwert von 21 % $_O2$ wird der Regelkreis deaktiviert, d.h. es wird auch keine Fehlerüberwachung durchgeführt. Das Gasmonitor-System (optional) ist trotzdem aktiv.

Hinweis

Bitte stellen Sie sicher, dass der Sensor korrekt in die Buchse eingesteckt wurde. Wenn der Sensor nicht korrekt installiert wurde, kann es zu Korrosion an den Kontakten und einer Fehlkalibrierung während der auto-start-Routine kommen. Sie können die Funktion einfach durch die Aktivierung des Sensors feststellen. Wenn nach 10 Minuten keine Fehlermeldung kommt, kann die auto-start-Routine durchgeführt werden.

auto-start-Funktion

Die Funktion auto-start ist eine automatisierte Routine zum Start und anschließenden Abgleich des $_{CO}$ 2-Messsystems. Nach dem Start regelt die Gerätesteuerung den eingestellten Temperatur-Sollwert ein. Gleichzeitig wird die Feuchte aufgebaut. Sind Temperatur und relative Feuchte auf einen konstanten Wert eingeregelt, wird das CO₂-Messsystem automatisch auf diese Werte abgeglichen und der Nutzraum mit dem voreingestellten CO₂-Wert begast.

Hinweise zur Anwendung der auto-start-Routine:

Damit die spezifizierte Genauigkeit des CO_2 -Messsystems erhalten bleibt, sollte das Gerät immer mit der auto-start-Routine gestartet werden, wenn:

- bei der Einstellung des Temperatur-Sollwertes ein Unterschied von mehr als 1 °C eingegeben wird,
- die Funktion Low-Feuchte aktiviert/deaktiviert wird,
- das Gerät nach längeren Betriebsunterbrechungen neu gestartet wird.

Im Rahmen der Reinigungs- und Wartungsarbeiten sollte die auto-start-Routine mindestens einmal vierteljährlich gefahren werden.

Dauer der auto-start-Routine:

Die auto-start-Routine dauert in der Regel 5 bis 7 Stunden. Bei niedrigen Raumtemperaturen und kaltem Gerät kann die auto-start-Routine bis zu 10 Std. dauern. Wird während der laufenden auto-start-Routine die Glastür geöffnet oder wird das Gerät vom Netz getrennt, wird die Routine unterbrochen und nach dem Schließen der Glastür oder dem Wiederanschluss an das Netz automatisch neu gestartet.

Startbedingungen für die auto-start-Routine:

Bei Beginn der auto-start-Routine darf die Atmosphäre im Nutzraum des Gerätes nur aus Umgebungsluft bestehen. Die Sollwerte für CO₂ und O₂ sollten vor dem Start auf die gewünschten Werte eingestellt werden. Das Wasserreservoir des Nutzraumes muss mit ausreichend Wasser befüllt werden.

Verhinderung des Starts der auto-start-Routine:

Die auto-start-Routine kann nicht gestartet werden, wenn einer der folgenden Fehler im Moment ansteht.

Regelkreis Temperatur:

- Sensor-/Fühlerbruch,
- Istwert über Sollwert (Abweichung zu stark),
- Istwert unter Sollwert (Abweichung zu stark),
- Istwert nicht plausibel
- Kalibrierwerte zu groß bzw. zu klein
- Keine Kommunikation zum Sensor

Regelkreis CO₂-Gasversorgung:

keine Kommunikation zum Sensor.
 Die auto-start -Taste wird in einem solchen Fehlerfall ausgeblendet und ist nicht verfügbar.

Fehlerhafter Abbruch der auto-start-Routine:

Die auto-start-Routine wird abgebrochen, wenn:

- ein Fehler im Temperatur-Regelkreis detektiert wird,
- ein Fehler im CO₂-Regelkreis detektiert wird,
- der Füllstand des Wassers zu niedrig ist,
- der einzustellende _{CO}2-Wert außerhalb der Toleranz liegt.

auto-start-Routine trocken ausführen:

Soll die auto-start-Routine trocken - ohne Wasserfüllung im Wasserreservoir des Nutzraumes ausgeführt werden, muss vor dem Start der Wasserstandsensor deaktiviert werden (Kapitel "Optionen" auf Seite 6-31).

auto-start aktivieren

Vorbereitungen zum Start:

- 1. Sicherstellen, dass die Ventile der CO₂-/O₂-/N₂-Gasversorgungsanlage geöffnet sind.
- 2. Ausreichend aufbereitetes Wasser in das Wasserreservoir des Nutzraumes gießen. Die Füllmenge darf die obere Füllstandsmarke nicht überschreiten.

auto-start-Routine aufrufen:

- 1. Taste auto-start drücken.
- Das Menü auto-start-Ablauf wird angezeigt.



Abbildung 6-13. auto-start aktivieren

- 2. Menü auto-start-Ablauf wieder verlassen und auto-start abbrechen:
- Taste Ende drücken.
- 3. Ggf. Wasser nachfüllen.
- 4. auto-start starten:
- Taste **Start** drücken.

- 5. Nutzraum belüften, beide Gerätetüren öffnen, bis nach 30 s das Zeitsignal ertönt.
- 6. Dann beide Gerätetüren schließen.
- Die Statusanzeige mit Daten zum Prozessverlauf wird angezeigt.



Abbildung 6-14. Statusanzeige auto-start-Status

Hinweis Abbruch: Die auto-start-Routine kann jederzeit unterbrochen werden! Taste **Stop** drücken.

Automatischer Neustart:

Die auto-start-Routine wird automatisch neu gestartet, wenn der Ablauf der Routine durch eine der folgenden Aktionen abgebrochen wurde:

- Öffnen der Glastür,
- Öffnen der Außentür bei optionaler Gasblende,
- Unterbrechung der Stromversorgung

auto-start unterbrechen

Wurde in der Statusanzeige die Taste **Stop** gedrückt, wird die auto-start-Routine unterbrochen und als Sicherheitsabfrage das Dialogfenster auto-start-Stop aufgerufen. Die Routine kann jetzt endgültig abgebrochen oder wieder fortgesetzt werden.

- 1. auto-start fortsetzen:
- Taste **Zurück** drücken.
- Es erfolgt ein Rücksprung in die Statusanzeige, die auto-start-Routine wird fortgesetzt.
- 2. auto-start abbrechen:
- Taste Ende drücken.
- Das Warndreieck wird als Abbruchhinweis zusammen mit dem akustischen Signal ausgegeben.



Abbildung 6-15. auto-start unterbrechen

- 3. Abbruchhinweis bestätigen:
- An eine beliebige Stelle des Displays drücken.
- Das Dialogfenster Fehler mit der Bezeichnung des Fehlers wird angezeigt.

Fehler		
System - F	ehler: auto-start	

Abbildung 6-16. Fehlermeldung nach Unterbrechen von auto-start

- 4. Fehlermeldung bestätigen:
- Taste **Ende** drücken.
- Es erfolgt ein Rücksprung ins Hauptmenü.

Steri-run aufrufen

Steri-run ist eine automatische Dekontaminationsroutine zur Desinfektion des Gerätenutzraumes. Die Steri-run-Dekontaminationsroutine dekontaminiert in einem automatisierten Programmablauf den kompletten Nutzraum inklusive des Regalsystems und der Sensoren. Ein detaillierter Ablauf der Gerätefunktion ist im Kapitel Reinigung und Desinfektion beschrieben (Kapitel 8, Seite 1).

Benutzerkonfiguration

Die Einstellungen zur Benutzerkonfiguration ermöglichen eine Anpassung der Benutzeroberfläche und der Gerätezusatzfunktionen an die Erfordernisse des alltäglichen Betriebes.Das Auswahlmenü zur Benutzerkonfiguration (Abbildung 6-17) wird über die Taste Menü des Hauptbildschirms aufgerufen.

Einstellung / Setup	Tastensperre
Datenaufzeichnung	Softwareversionen
Optionen	1
lconerklärung	1

Abbildung 6-17. Menü Benutzerkonfiguration

Das Menü Benutzerkonfiguration ist in sechs Kategorien gegliedert:

- Einstellungen/Setup,
- Datenaufzeichnung,
- Optionen,
- Iconerklärung,
- Tastensperre,
- Softwareversionen.

Um eine benutzerspezifische Einstellung in einem Dialogfenster vorzunehmen, wird durch die in den Abbildungen angegebenen Untermenüs navigiert und dann das Dialogfenster aufgerufen.

Einstellungen/Setup

Das Auswahlmenü **Einstellungen/Setup** (Abbildung 6-18) ermöglicht den Zugriff auf eine Reihe von Eingabedialogen mit Einstellmöglichkeiten zur individuellen Konfiguration der Benutzeroberfläche und der Benutzerschnittstelle:

- Code für Tastensperre ändern,
- Datum / Uhrzeit einstellen,
- Helligkeit des Displays einstellen,
- Tastenton einstellen,
- Schnittstelle einrichten,
- Sprache Bedienoberfläche einstellen,
- Erinnerungsintervalle einstellen.

Tastensperre Code	Schnittstelle
Datum/Uhrzeit	Sprache
Helligkeit	Erinnerungsintervalle
Tastenton	ב
	Ende

Abbildung 6-18. Menü Einstellungen/Setup

Code für Tastensperre ändern

Die Tastensperre verhindert eine unbefugte Änderung der Betriebseinstellungen. Es werden dadurch nur solche Tasten gesperrt, die eine Werteingabe bewirken.

Der Code der Tastensperre besteht aus einer Pflichteingabe von vier Ziffern.

• Die werksseitige Vorgabe ist: 0000.

Diese Vorgabe kann auf einen benutzerdefinierten Code geändert werden und anschließend mit dem Dialogfenster TASTENSPERRE ("Tastensperre aktivieren/deaktivieren" auf Seite 6-42) aktiviert werden.

Code für Tastensperre ändern:

- 1. Taste Menü drücken.
- 2. Menübefehl Tastensperre Code wählen.
- Es erscheint der in Abbildung 6-19 dargestellte Eingabedialog.

Tastensperre Code	Code
	6 7 8 9 0
	Löschen Enter Zurück

Abbildung 6-19. Code für Tastensperre ändern

- 3. Werksvorgabe 0000 eingeben:
- Entsprechende Zifferntasten drücken.
- Die Ziffernkombination wird im Eingabefeld verdeckt angezeigt.

- 4. Eingabe bestätigen:
- Taste Enter drücken.
- In der Anzeige erscheint die Eingabeaufforderung Code Neu. Vier Leerstellen zeigen an, dass der neue Tastencode eingegeben werden kann.

Code neu		
1 2 3 4 6 7 8 9	5	
Enter	<<	

Abbildung 6-20. Code für Tastensperre ändern

Neuen 4-stelligen Code eingeben:

- Entsprechende Zifferntasten drücken.
- Die Ziffernkombination wird im Eingabefeld angezeigt.
- 5. Cursor zum Überschreiben eines Wertes nach links setzen:
- Taste **Rückschritt** (<<) drücken.
- 6. Eingabe übernehmen und speichern:
- Taste Enter drücken.
- Es erfolgt ein Rücksprung ins Auswahlmenü Einstellungen/Setup.
- In der Icon-Leiste im Hauptmenü (Abbildung 6-2 auf Seite 3) erscheint das Schlosssymbol als Aktivitätsanzeige für die Tastensperre.

Hinweis Benutzerdefinierten Code ändern:

Der benutzerdefinierte Code kann mit derselben Verfahrensweise beliebig oft verändert werden:

- Neucodierung durch Eingabe des gültigen Codes aktivieren,
- neuen Code eingeben und bestätigen.

Datum/Uhrzeit einstellen

Der Eingabe-Dialog ermöglicht die Einstellung von Datum und Uhrzeit auf die gewünschte Zeitzone.

- 1. Taste **Menü** drücken.
- 2. Menübefehl Datum / Uhrzeit wählen.
- Es erscheint der in Abbildung 6-21 dargestellte Auswahldialog.

Datum/Uhrzeit	Datum
	Ende

Abbildung 6-21. Auswahlmenü Datum/Uhrzeit

- 3. Zum Ändern des Datums die Option Datum wählen.
- Es erscheint der in Abbildung 6-22 dargestellte Eingabedialog.

Datum			
	Datum	05.10.13	
		3 4 5	
	6 7	8 9 0	
		nter <<)

Abbildung 6-22. Datum einstellen

- 4. Datum eingeben:
- Zifferntasten drücken.
- Die eingegebenen Ziffern werden im Eingabefeld angezeigt.
- 5. Cursor zum Überschreiben eines Wertes nach links setzen:
- Taste **Rückschritt** (<<) drücken.
- 6. Eingabe übernehmen und speichern:
- Taste Enter drücken.
- 7. Es erfolgt ein Rücksprung ins Auswahlmenü Datum / Uhrzeit.
- 8. Zum Ändern der Uhrzeit die Option Uhrzeit wählen.
- Es erscheint der in Abbildung 6-23 dargestellte Eingabedialog.



Abbildung 6-23. Uhrzeit einstellen

- 9. Uhrzeit eingeben:
- Zifferntasten drücken.
- Die eingegebenen Ziffern werden im Eingabefeld angezeigt.
- 10. Cursor zum Überschreiben eines Wertes nach links setzen:
 - Taste **Rückschritt** (<<) drücken.
- 11. Eingabe übernehmen und speichern:
- Taste Enter drücken.

12. Es erfolgt ein Rücksprung ins Auswahlmenü Datum / Uhrzeit.

Helligkeit des Displays einstellen

Der Eingabedialog ermöglicht die Einstellung der Helligkeit des Bedienfeldes im Wertebereich von 1 bis 100 %.

Helligkeit einstellen

- 1. Taste Menü drücken.
- 2. Menübefehl Einstellungen / Setup wählen.
- 3. Option Display Helligkeit wählen.
- Es erscheint der in Abbildung 6-24 dargestellte Eingabedialog.

Helligkeit	Display Hell	igkeit		
	aktuell	31		
	neu	31	%	-
	·	Enter	+	כ

Abbildung 6-24. Displayhelligkeit einstellen

- 4. Wert erhöhen:
- Taste + drücken.
- 5. Wert verringern:
- Taste drücken.
- 6. Die Wertänderung wird im Anzeigefeld angezeigt. Der Hinweis **neu** zeigt an, dass der Wert geändert, jedoch noch nicht gespeichert wurde.
- 7. Änderung übernehmen und speichern:
- Taste Enter drücken.
- Es erfolgt ein Rücksprung ins Auswahlmenü Einstellungen/Setup.

Tastenton einstellen

Der Eingabedialog ermöglicht die Einstellung der Lautstärke des Tastentones, der beim Betätigen einer Taste ausgegeben wird.

Der Wertebereich reicht von 0 bis 100. Die Veränderung erfolgt in 5 %-Schritten.

Lautstärke des Tastentons einstellen

- 1. Taste Menü drücken.
- 2. Menübefehl Einstellungen / Setup wählen.
- 3. Option Tastenton wählen.
- Es erscheint der in Abbildung 6-25 dargestellte Eingabedialog.
| Tastenton | Lautstärke T | astenton | | D |
|-----------|--------------|----------|---|---|
| | aktuell | ₅
5 | % | |
| | · · | Enter | + | |

Abbildung 6-25. Lautstärke des Tastentons einstellen

- 1. Wert erhöhen:
- Taste + drücken.
- 2. Wert verringern:
- Taste drücken.
- 3. Die Wertänderung wird im Anzeigefeld angezeigt. Der Hinweis **neu** zeigt an, dass der Wert geändert, jedoch noch nicht gespeichert wurde.
- 4. Änderung übernehmen und speichern:
- Taste Enter drücken.
- Es erfolgt ein Rücksprung ins Auswahlmenü Einstellungen/Setup.

Baudrate USB-Schnittstelle einstellen

Der Eingabedialog ermöglicht die Einstellung der Datenübertragungsrate der USB-Schnittstelle:

Die Datenübertragungsrate der Schnittstelle kann innerhalb der definierten Baudraten (9600, 19200, 38400, 57600 Baud) verändert werden.

Baudrate einstellen

- 1. Taste Menü drücken.
- 2. Menübefehl Einstellungen / Setup wählen.
- Es erscheint der in Abbildung 6-26 dargestellte Auswahldialog.

Schnittstelle	Schnittstelle USB
	Ende

Abbildung 6-26. Auswahlmenü Schnittstelle USB

- 3. Option Schnittstelle USB wählen (Abbildung 6-26).
- Es erscheint der in Abbildung 6-27 dargestellte Eingabedialog.

Schnittstelle USB	Baudrate	USB	
	aktuell	57600	
	neu	57600	Baud
	_ ·	Enter	+

Abbildung 6-27. Baudrate USB-Schnittstelle einstellen

- 4. Mit Taste + oder Baudrate der USB-Schnittstelle anpassen (Abbildung 6-27).
- Baudrate erhöhen: Taste + drücken.
- Baudrate verringern: Taste drücken.
- 5. Die Wertänderung wird im Anzeigefeld angezeigt. Der Hinweis **neu** zeigt an, dass der Wert geändert, jedoch noch nicht gespeichert wurde.
- 6. Änderung übernehmen und speichern:
- Taste Enter drücken.
- Es erfolgt ein Rücksprung ins Auswahlmenü Schnittstelle USB.
- 7. Neue Einstellungen aktivieren:
- Zurückblättern ins Hauptmenü.
- Ca. 10 s warten und einen Neustart durchführen, Gerät am Netzschalter aus-/einschalten.

Sprache der Bedienoberfläche einstellen

Der Eingabedialog ermöglicht die Einstellung der Sprache der Bedienoberfläche. Es stehen sieben Sprachvarianten zur Auswahl:

- deutsch,
- englisch,
- spanisch,
- französisch,
- italienisch,
- chinesisch,
- japanisch.

Anzeigesprache auswählen:

- 1. Taste Menü drücken.
- 2. Menübefehl Sprache wählen.
- Es erscheint der in Abbildung 6-28 dargestellte Eingabedialog.

Sprache			
	 Sprachei	instellung	
	aktuell	deutsch	
	neu	deutsch	
	-	Enter	+

Abbildung 6-28. Sprache einstellen

- 3. In der Auswahl nach oben blättern:
- Taste + drücken.
- 4. In der Auswahl nach unten blättern:
- Taste drücken.
- 5. Die neu gewählte Sprache wird im Anzeigefeld angezeigt. Der Hinweis **neu** zeigt an, dass der Wert geändert, jedoch noch nicht gespeichert wurde.
- 6. Auswahl übernehmen und speichern:
- Taste Enter drücken.
- Es erfolgt ein Rücksprung ins Auswahlmenü Einstellungen/Setup.

Erinnerungsintervalle einstellen

Die Erinnerungsintervalle sind Teil des Melde- und Kontrollsystems der Gerätesteuerung. Für die beiden wichtigen Gerätefunktionen Steri-run und auto-start, sowie für routinemäßige Servicearbeiten können jeweils Zeitpunkte bestimmt werden, bei deren Fälligkeit eine Meldung ausgelöst wird. Zählbeginn ist 00:00 Uhr des Tages, an dem das jeweils zuvor eingestellte Erinnerungsintervall abgelaufen ist.

Am Tage der Fälligkeit wird für das aktivierte Erinnerungsintervall eine Hinweismeldung am Display ausgegeben:

- Steri-run: Bitte Steri-run ausführen.
- auto-start: Bitte auto-start ausführen. Erscheint nach erfolgreich durchlaufener Steri-run-Dekontaminationsroutine.
- Serviceintervall: Bitte Service durchführen. Die Servicemeldung kann bestätigt werden. In der Folge wird das Icon Service anfordern angezeigt.

Die Hinweismeldungen werden ausgeblendet, wenn die Routinen erfolgreich durchgeführt wurden.

Werkseitige Einstellungen

Steri-run-Dekontaminationsroutine	90 Tage
auto-start-Routine	Aus
Serviceintervall	Aus
HEPA-Filter-Intervall	365 Tage

Erinnerungsintervalle einstellen

- 1. Taste Menü drücken.
- 2. Menübefehl Erinnerungsintervalle wählen.
- Es erscheint der in Abbildung 6-29 dargestellte Auswahldialog.

Erinnerungsintervalle	steri-run Rücksetzen HEPA	
	auto-start	-
	Serviceintervall	
	HEPA-Filter-Intervall	
	Ende]

Abbildung 6-29. Funktion für Erinnerungsintervall auswählen

- 3. Entsprechenden Menübefehl wählen, z.B. steri-run.
- Es erscheint der in Abbildung 6-29 dargestellte Eingabedialog.

steri-run	Erinnerungs	sintervall steri-run	
	aktuell	365	
	neu	100	Tage
	· ·	Enter	+

Abbildung 6-30. Erinnerungsintervall für steri-run einstellen

- 1. Anzahl der Tage erhöhen:
- Taste + drücken.
- 2. Anzahl der Tage verringern:
- Taste drücken.
- 3. Die Wertänderung wird im Anzeigefeld angezeigt. Der Hinweis **neu** zeigt an, dass der Wert geändert, jedoch noch nicht gespeichert wurde.
- 4. Erinnerungsintervall deaktivieren:
- Wert auf **Aus** setzen.
- Taste drücken.
- 5. Änderung übernehmen und speichern:
- Taste Enter drücken.
- Es erfolgt ein Rücksprung ins Auswahlmenü Erinnerungsintervalle.

Datenaufzeichnung

Das Auswahlmenü Datenaufzeichnung (Abbildung 6-31) ermöglicht den Zugriff auf Dialoge zur Aufzeichnung und Ausgabe von Ereignissen während des Gerätebetriebes:

- Ereignisanzeige,
- Zeitintervall (des Speicherzyklus),
- Fehlertabelle.

Datenaufzeichnung	
	Speicherzykluszeit
	Fehlertabelle
	Ende

Abbildung 6-31. Auswahlmenü Datenaufzeichnung

Ereignisse anzeigen

Die Ereignisanzeige berichtet in Form von kurzen einzeiligen Einträgen unter Angabe von Datum und Uhrzeit über die aufgezeichneten Ereignisse im Gerätebetrieb.

Die Einträge werden in zeitlicher Reihenfolge - mit dem jüngsten Ereignis an oberster Position gelistet. Die Liste kann angesehen, aber nicht editiert werden. Sofern die Ereignisanzeige aus mehreren Seiten besteht, kann in der Liste geblättert werden. Die Statusanzeige gibt Auskunft darüber, welche Seite der Gesamtseitenzahl aktuell angezeigt wird.

Ereignisanzeige aufrufen

- 1. Taste Menü drücken.
- 2. Menübefehl Datenaufzeichnung wählen.
- Es erscheint der in Abbildung 6-31 dargestellte Auswahldialog.
- 3. Menübefehl Ereignisanzeige wählen.
- Es erscheint die in Abbildung 6-32 dargestellte Listenanzeige.





- 1. In der Liste vorblättern:
- Taste **Weiter** drücken.
- 2. In der Liste zurückblättern:
- Taste Vorher drücken.
- 3. Anzeige beenden:
- Taste Ende drücken.
- Es erfolgt ein Rücksprung ins Auswahlmenü Datenaufzeichnung.

Speicherzyklus einstellen:

Auf Grund von limitierten Speicherressourcen werden beim Aufzeichnen von neuen Ereignissen die ältesten Einträge gelöscht. Der Zeitraum, aus dem die angezeigten Ereignisse stammen, hängt maßgeblich von der gewählten Speicherzykluszeit ab.

Ereignisanzeige aufrufen

- 1. Taste Menü drücken.
- 2. Menübefehl Datenaufzeichnung wählen.
- Es erscheint der in Abbildung 6-31 dargestellte Auswahldialog.
- 3. Menübefehl Speicherzykluszeit wählen.
- Es erscheint der in Abbildung 6-33 dargestellte Eingabedialog.



Abbildung 6-33. Speicherzyklus einstellen

Speicherzyklus	Maximal angezeigter Zeitraum
10 s	22,5 Stunden
30 s	2,8 Tage
60 s	5,6 Tage
120 s	11,2 Tage
180 s	16,8 Tage
3600 s	336 Tage

Die Einstellung steuert den Speicherzyklus in Sekundenabschnitten, in denen die Messwerte der Regelkreise im Gerätebetrieb aufgezeichnet und in der Verlaufsanzeige wiedergegeben werden ("Verlaufsanzeige skalieren" auf Seite 6-43).

Die Einstellungen können im Wertebereich von 10 s bis 3600 s vorgenommen werden.

- 1. Wert erhöhen:
- Taste + drücken.
- 2. Wert verringern:
- Taste drücken.
- 3. Die Wertänderung wird im Anzeigefeld angezeigt. Der Hinweis **neu** zeigt an, dass der Wert geändert, jedoch noch nicht gespeichert wurde.
- 4. Änderung übernehmen und speichern:
- Taste Enter drücken.
- Es erfolgt ein Rücksprung ins Auswahlmenü Datenaufzeichnung.

Hinweis Speicherzyklen der Datenaufzeichnung: Die Speicherzykluszeit hat keinen Einfluss auf die Einträge der Fehlertabelle.

Fehlertabelle einsehen

Die Fehlertabelle listet die vom geräteinternen Kontrollsystem detektierten Fehler in absteigender zeitlicher Reihenfolge. Der zuletzt detektierte Fehler wird an die oberste Position von insgesamt 22 möglichen Einträgen gesetzt. Ein Eintrag besteht aus dem Regelkreis, in dem der Fehler detektiert wurde, Datum, Uhrzeit und einer Fehlerbezeichnung. Die Fehlertabelle kann eingesehen, aber nicht editiert werden. Sofern die Ereignisanzeige aus zwei Seiten besteht, kann in der Liste geblättert werden. Die Statusanzeige **001/002 bzw. 002/002** gibt Auskunft darüber, welche der beiden Seiten aktuell angezeigt wird.

Fehlertabelle einsehen

- 1. Taste Menü drücken.
- 2. Menübefehl Datenaufzeichnung wählen.
- Es erscheint der in Abbildung 6-31 dargestellte Auswahldialog.
- 3. Menübefehl Ereignisanzeige wählen.
- Es erscheint die in Abbildung 6-34 dargestellte Listenanzeige.

Fehlertabelle				
	Rkreis	Datum	Uhrzeit	Fehler 001/002
	CO2	05.10.13	20:59:51	Störung Kommunikation
	RH	05.10.13	00:21:46	kein Wasser
	CO2	28.09.13	00:31:32	Störung Kommunikation
	RH	28.09.13	00:31:19	kein Wasser
	SYS	16.08.13	11:31:39	IR-Sensor getauscht
	SYS	16.08.13	11:30:47	Fehler EEPROM Mainboard
	RH	16.08.13	11:12:13	kein Wasser
	Т	16.08.13	11:12:11	Sensor-/Fühlerbruch
	RH	16.08.13	11:11:59	kein Wasser
	RH	16.08.13	11:10:08	kein Wasser
	т	16.08.13	11:09:56	Sensor-/Fühlerbruch
			End	le Weiter

Abbildung 6-34. Fehlertabelle einsehen

Hinweis Fehlerbehandlung:

Eine ausführliche Übersicht zu Fehlerursachen und Fehlerbeseitigung finden Sie am Ende dieses Kapitels!

- 1. In der Fehlertabelle vorblättern:
- Taste Weiter drücken.
- 2. In der Liste zurückblättern:
- Taste Vorher drücken.
- 3. Anzeige beenden:
- Taste **Ende** drücken.
- Es erfolgt ein Rücksprung ins Auswahlmenü Datenaufzeichnung.

Optionen

Das Auswahlmenü **Optionen** (Abbildung 6-35) ermöglicht den Zugriff auf alle Einstellungsdialoge für die funktionalen Geräteoptionen:

- Alarm,
- Low Feuchte,
- Gasblende (optional),
- Luftfeuchte-Sensor,
- Hupe,
- O₂-Gasversorgung (optional),
- HEPA-Filter.

Optionen		
	Alarm	Hupe
	Low Feuchte	02
	Gasblende	HEPA-Filter
	Wasserstandsensor	
		inde



Alarmrelais einstellen

Das Alarmrelais bildet die Schnittstelle für die Anbindung des geräteinternen Kontrollsystems an ein externes Monitoring-System zur Überwachung der Stromversorgung. Abhängig vom erforderlichen Eingangssignal des externen Monitoring-Systems kann die Netzüberwachung ein- oder ausgeschaltet werden. Ist die Netzüberwachung eingeschaltet, wird ein Stromausfall als Fehler erkannt. Die Verdrahtung des Alarmrelais wird im Abschnitt "Alarmkontakt anschließen:" auf Seite 4-18 beschrieben.

- 1. Taste Menü drücken.
- 2. Menübefehl **Optionen** wählen.
- Es erscheint der in Abbildung 6-35 dargestellte Auswahldialog.
- 3. Menübefehl Alarm wählen.
- Es erscheint der in Abbildung 6-36 dargestellte Auswahldialog.

Alarm	Alarmrelais
	Ende

Abbildung 6-36. Auswahlmenü Alarm

- 4. Menübefehl Alarmrelais wählen.
- Es erscheint der in Abbildung 6-37 dargestellte Eingabedialog.

Alarmrelais	Alarmrelais - Netzüberwachung
	aktueli an
	neu an
	- Enter +

Abbildung 6-37. Alarmrelais einstellen

- 1. Zwischen zwei Zuständen hin- und herschalten:
- Taste + drücken.
 - oder
- Taste drücken.
- 2. Änderung übernehmen und speichern:
- Taste Enter drücken.
- Es erfolgt ein Rücksprung ins Auswahlmenü Optionen.

Low Feuchte einstellen

Bildet sich aufgrund der hohen relativen Feuchte an den Kulturgefäßen ein Tauniederschlag, kann die Feuchte im Nutzraum auf einen niedrigeren Wert angepasst werden. Werkseitig ist die

Gerätessteuerung auf High-Feuchte - ca. 93 % relative Feuchte- voreingestellt.

Mit der Aktivierung der Low Feuchte wird die relative Feuchte im Nutzraum von ca. 93 % auf ca. 90 % abgesenkt.

Die Veränderung beansprucht eine lange Anpassungsphase. Damit sie wirksam eine Betauung von Kulturgefäßen verhindert, muss sie als Dauereinstellung angewendet werden.

Luftfeuchte im Nutzraum senken

- 1. Taste Menü drücken.
- 2. Menübefehl Optionen wählen.
- Es erscheint der in Abbildung 6-35 dargestellte Auswahldialog.
- 3. Menübefehl Low Feuchte wählen.
- Es erscheint der in Abbildung 6-38 dargestellte Eingabedialog.

Low Feuchte			
	Low Feucht	e	
	aktuell	aus	
	neu	aus	
	-	Enter	+

Abbildung 6-38. Low Feuchte einstellen

- 1. Zwischen zwei Zuständen hin- und herschalten:
- Taste + drücken. oder
- Taste drücken.
- 2. Änderung übernehmen und speichern:
- Taste Enter drücken.

Es erfolgt ein Rücksprung ins Auswahlmenü Optionen.



Nach der Rückkehr zum Hauptmenü erscheint das Icon Low Feuchte.

Hinweis Low-Feuchte:

Das Ein-/Ausschalten der Funktion Low-Feuchte wird in die Ereignisliste eingetragen.

Gasblende einstellen

Weil die Öffnungsquerschnitte beim Zugriff auf die Proben kleiner sind, erreichen Geräte, die mit der optionalen Gasblende ausgerüstet sind, kürzere Erholungszeiten der Inkubationsparameter:

- Nutzraumtemperatur,
- CO₂-Konzentration,
- O₂-Konzentration,
- Relative Feuchte.

Nach einer Umrüstung des Gerätes muss die Gerätesteuerung auf die Option Gasblende umgestellt werden.

Hinweis Fehlfunktion:

Die Umstellung auf Gasblende bewirkt eine Änderung der Steuerungsparameter. Wird die Funktion Gasblende nicht entsprechend der tatsächlich eingebauten Türvariante eingestellt, kann es zu Fehlfunktionen bei der Inkubationsleistung kommen.

Gasblende einstellen

- 1. Taste **Menü** drücken.
- 2. Menübefehl **Optionen** wählen.
- Es erscheint der in Abbildung 6-35 dargestellte Auswahldialog.

- 3. Menübefehl Gasblende wählen.
- Es erscheint der in Abbildung 6-39 dargestellte Eingabedialog.

Gasblende			
	Gasblende		
	aktuell	Glastür	
	neu	Glastür	
	-	Enter	+

Abbildung 6-39. Gasblende einstellen

- 1. Zwischen zwei Optionen hin- und herschalten:
- Taste + drücken.
- oder
- Taste drücken.
- 2. Änderung übernehmen und speichern:
- Taste Enter drücken.
- Es erfolgt ein Rücksprung ins Auswahlmenü Optionen.

Wasserstandsensor ein- / ausschalten

Für einen Inkubationsbetrieb bei Umgebungsfeuchte oder, wenn die auto-start-Routine trocken - ohne Wasserfüllung - ausgeführt werden soll, kann der Wasserstand-Sensor ausgeschaltet werden. Damit werden Alarmmeldungen des Wasserstandsensors durch das geräteinterne Kontrollsystem unterbunden:



WARNUNG Wenn der Wasserstandsensor ausgeschaltet ist, kann die Steri-run-Routine trotz Wasser im Wasserreservoir gestartet werden. Dies stellt eine Fehlbenutzung dar und kann den Lüftermotor zerstören.

- 1. Taste Menü drücken.
- 2. Menübefehl Optionen wählen.
- Es erscheint der in Abbildung 6-35 dargestellte Auswahldialog.
- 3. Menübefehl Wasserstandsensor wählen.
- Es erscheint der in Abbildung 6-40 dargestellte Eingabedialog.

Wasserstandsensor			
	Wasserstan	dsensor	
	aktuell	an	
	neu	an	
	-	Enter	+

Abbildung 6-40. Wasserstandsensor einstellen

- 1. Zwischen zwei Zuständen hin- und herschalten:
- Taste + drücken. oder
- Taste drücken.
- 2. Änderung übernehmen und speichern:
- Taste Enter drücken.
- Es erfolgt ein Rücksprung ins Auswahlmenü Optionen.

Hupe ein- / ausschalten:

Hat das geräteinterne Kontrollsystem einen Fehler erkannt, wird:

- zusätzlich zur optischen Fehlermeldung und zur Schaltung des Alarmrelais,
- ein akustischer Alarm als abgesetzter Hupton ausgegeben.

Der Hupton kann dauerhaft ausgeschaltet werden.

- 1. Taste **Menü** drücken.
- 2. Menübefehl **Optionen** wählen.
- Es erscheint der in Abbildung 6-35 dargestellte Auswahldialog.
- 3. Menübefehl Hupe wählen.
- Es erscheint der in Abbildung 6-41 dargestellte Eingabedialog.

Нире				
		Akustischer	Alarm	
		aktuell	an	
		neu	an	
	(_	Enter	•

Abbildung 6-41. Alarmrelais einstellen

- 1. Zwischen zwei Zuständen hin- und herschalten:
- Taste + drücken.
 - oder
- Taste drücken.
- 2. Änderung übernehmen und speichern:
- Taste Enter drücken.
- Es erfolgt ein Rücksprung ins Auswahlmenü Optionen.

O₂-Regelung ein-/ausschalten

Abhängig von den Anforderungen an den Arbeitsprozess kann die O₂-Regelung ein- bzw. ausgeschaltet werden. Diese Einstellung ist nur bei Ausstattung mit optionaler O₂-/N₂-Regelung möglich.

- 1. Taste **Menü** drücken.
- 2. Menübefehl **Optionen** wählen.
- Es erscheint der in Abbildung 6-35 dargestellte Auswahldialog.
- 3. Menübefehl O₂ wählen.
- Es erscheint der in Abbildung 6-42 dargestellte Eingabedialog.

O2-Konfigur	ation	
aktuell	an	
neu	an	
·	Enter	+

Abbildung 6-42. O₂-Regelung ein-/ausschalten

- 1. Zwischen zwei Zuständen der O2-Regelung hin- und herschalten:
- Taste + drücken. oder
- Taste drücken.
- 2. Die Wertänderung wird im Anzeigefeld angezeigt. Der Hinweis neu zeigt an, dass der Wert geändert, jedoch noch nicht gespeichert wurde.
- 3. Einstellung übernehmen und speichern:
- Taste Enter drücken.
- Es erfolgt ein Rücksprung ins Auswahlmenü Optionen.

Hinweis Anzeige O₂-Wert:

Ist die O₂-Regelung ausgeschaltet, wird im O₂-Anzeigefeld kein Istwert (- - -) angezeigt. Diese Verfahrensweise wirkt sich schonend auf die Beanspruchung des O₂-Sensors aus. Ist der Sollwert auf 21 % eingestellt, findet keine Überwachung des O₂-Regelkreises statt. Dies gilt für beide O₂-Regelbereiche:

- Regelbereich I: 1 % 21 %
- Regelbereich II: 5 % 90 %

Im O2-Anzeigefeld wird dann der Istwert angezeigt.

Nutzraum lüften

Wurde das Gerät mit O_2 oder N_2 betrieben, muss nach dem Ausschalten der O_2 -Regelung der Nutzraum gelüftet werden.

Gasmonitoring

Der Schaltzustand des O₂-Regelkreises hat keinen Einfluss auf die Funktion des optionalen Gasmonitorings. Auch bei ausgeschalteter O₂-Regelung oder deaktiviertem O₂-Sensor ist das Gasmonitor-System aktiv.

HEPA-Filter aktivieren / deaktivieren:

Wenn das Gerät ohne den eingebauten HEPA-Filter betrieben werden soll, muss dieser in der Konfiguration zur Vermeidung von Fehlfunktionen deaktiviert werden.

- 1. Taste Menü drücken.
- 2. Menübefehl Optionen wählen.
- Es erscheint der in Abbildung 6-35 dargestellte Auswahldialog.
- 3. Menübefehl HEPA-Filter wählen.
- Es erscheint der in Abbildung 6-43 dargestellte Auswahldialog.



Abbildung 6-43. HEPA-Konfiguration

- 4. Menübefehl HEPA-Konfiguration wählen.
- Es erscheint der in Abbildung 6-44 dargestellte Eingabedialog.

HEPA-Konfiguration	HEPA-Filter		
	aktuell	an	
	neu	an	
	-	Enter	+

Abbildung 6-44. HEPA-Filter aktivieren / deaktivieren

- 1. Zwischen zwei Zuständen hin- und herschalten:
- Taste + drücken. oder
- Taste drücken.

- 2. Änderung übernehmen und speichern:
- Taste Enter drücken.
- Es erfolgt ein Rücksprung ins Auswahlmenü Optionen.
- In der Icon-Leiste im Hauptmenü (Abbildung 6-2 auf Seite 3) erscheint nach 5 Minuten die Aktivitätsanzeige **ISO 5** für den HEPA-Filter.

Iconerklärung

ISO

5

Wichtige Betriebszustände oder Fehlermeldungen, z.B. die Tastensperre oder die Low Feuchte, werden zusätzlich zu den Eintragungen in der Datenaufzeichnung oder in der Fehlertabelle, als Icon im Hauptmenü des Tastbildschirms angezeigt. Die Bedeutung der Icons ist im Dialogfenster Iconerklärung erklärt.

Iconerklärung aufrufen

- 1. Taste Menü drücken.
- 2. Menübefehl Iconerklärung wählen.
- Es erscheint der in Abbildung 6-45 dargestellte Informationsdialog.



Abbildung 6-45. lconerklärung

- Anzeige beenden:
- Taste **Ende** drücken.
- Es erfolgt ein Rücksprung ins Auswahlmenü Benutzerkonfiguration.

Die funktionale Bedeutung der einzelnen Icons:

Gasflasche leer:



Fehleranzeige, die darauf hinweist, dass der Füllstand einer oder mehrerer Gasflaschen zu niedrig ist, um eine reguläre Gasversorgung zu gewährleisten.

Die Überwachungsfunktion ist nur bei Ausstattung des Gerätes mit einem Gasmonitoring-System (optional) verfügbar.

Übertemperatur:



Fehleranzeige, die darauf hinweist, dass die Gerätesteuerung den Übertemperaturschutz aktiviert hat und auf eine Notregelung umgeschaltet hat.

Tastensperre:



Funktionsanzeige, die anzeigt, dass die Tastensperre aktiviert wurde und aktuell keine Änderungen an den Einstellungen möglich sind (Anleitung siehe "Tastensperre aktivieren/deaktivieren" auf Seite 6-42).

Low Feuchte:



Funktionsanzeige, die anzeigt, dass die relative Feuchte im Nutzraum von ca. 93 % auf ca. 90 % abgesenkt wurde (Anleitung siehe "Low Feuchte einstellen" auf Seite 6-33).

HEPA-Filter aktiv:



Funktionsanzeige, die anzeigt, dass das HEPA-Filter im Nutzraum aktiviert wurde (Anleitung siehe "HEPA-Filter aktivieren / deaktivieren:" auf Seite 6-39).

Service anfordern:



Hinweis, der anzeigt, dass das routinemäßige Serviceintervall fällig ist. Die Anzeige des Icons wird durch die Zeitangabe im Dialogfenster Erinnerungsintervall gesteuert und wird angezeigt, nachdem die Erinnerungsmeldung bestätigt wurde.

Anzeige Füllstand Gasflaschen (optional):

Ist das Gerät mit der Funktion Gasmonitoring (optional) ausgestattet, werden die Icons für Gasflasche A und Gasflasche B im jeweiligen CO_2 - bzw. O_2 -Menü eingeblendet. Die Icons zeigen den Füllstand (voll/leer) der Gasflaschen an. Das blau gerahmte Icon zeigt dabei an, auf welche Gasflasche zur Weiterversorgung umgeschaltet werden kann.

Gasflaschenmo	nitor	Verlauf anzeigen
C02	4.3	0.0
Kalibration	Ende	Sollwert

Abbildung 6-46. Icons Gasmonitoring

- Der Wechsel auf eine volle Gasflasche kann manuell ausgeführt werden.
- Blau gerahmtes Flaschen-Icon drücken oder
- wird automatisch ausgeführt, wenn:
- der Flaschendruck unter 0,6 bar fällt.

Nach einem manuellen oder automatischen Wechsel der Gaszufuhr kann 30 s lang keine weitere Umschaltung vorgenommen werden. Ungefähr weitere 2 min dauert es, bis der Füllstand der Wechselflasche neu ermittelt ist und angezeigt wird.

Das Gasmonitoring-System überwacht den Füllstand beider angeschlossenen Gasflaschen. Ist eine Flasche leer wird:

- kein akustischer Alarm ausgegeben,
- eine Fehlermeldung angezeigt,
- ein Eintrag in die Ereignisliste geschrieben.

Sind beide Flaschen leer, wird:

- ein akustischer Alarm ausgegeben und das Alarmrelais wird geschaltet,
- eine Fehlermeldung angezeigt,
- ein Eintrag in die Fehlertabelle geschrieben,
- ein Eintrag in die Ereignisliste geschrieben.

Hinweis Flaschenwechsel:

Das automatische oder manuelle Umschalten zwischen den Flaschen wird in die Ereignisliste eingetragen.

Tastensperre aktivieren/deaktivieren

Der Eingabe-Dialog ermöglicht die Aufhebung bzw. die Aktivierung der Tastensperre. Werkseitig ist die Tastensperre auf den Standard-Code 0000 eingestellt.

- 1. Vierstelligen Code über die Zifferntasten eingeben. Die Eingabe wird im Anzeigefeld verschlüsselt angezeigt.
- 2. Falsche Eingabe komplett löschen:
- Taste Löschen drücken.
- 3. Eingabe abbrechen:
- Taste **Zurück** drücken.
- Es erfolgt ein Rücksprung ins Auswahlmenü Benutzerkonfiguration.
- 4. Eingabe bestätigen:
- Taste **Ende** drücken.
- Es erfolgt ein Rücksprung ins Auswahlmenü Benutzerkonfiguration.

Hinweis Bestehenden Code ändern:

Der aktuell gültige Code kann im Dialogfenster Tastensperre/Code des Menüs Einstellungen / Setup neu definiert werden ("Code für Tastensperre ändern" auf Seite 6-18).

Rücksetzung Code:

Ist der Code der Tastensperre nicht mehr verfügbar, kann die Rücksetzung auf den Standard-Code nur durch den Technischen Support von Thermo Fisher Scientific erfolgen.

Tastensperre		
	Code:	
		-
	Löschen Enter Zurück)

Abbildung 6-47. Tastensperre aktivieren/deaktivieren

Softwareversionen

Zeigt im Anzeigefeld die zur Gerätesteuerung implementierten Versionen der Software an.

- Anzeige beenden:
- Taste Ende drücken.
- Es erfolgt ein Rücksprung ins Auswahlmenü Benutzerkonfiguration.



Abbildung 6-48. Softwareversionen

Verlaufsanzeige skalieren

Die Verlaufsanzeige der drei Regelkreise:

- Temperatur,
- 0...20% CO₂,
- 0...100% O₂,

kann auf zwei unterschiedliche Darstellungsweisen skaliert werden.

- 1. Vollbild-Anzeige
- Taste **CO₂-Anzeigefeld** im Hauptmenü drücken.

• Das CO₂-Menü (Abbildung 6-9) wird angezeigt.



Abbildung 6-49. Verlaufsanzeige für CO₂-Konzentration aufrufen

- 2. Icon Verlauf anzeigen drücken.
- Die Verlaufsanzeige wird eingeblendet.



Abbildung 6-50. Verlaufsanzeige für CO₂-Konzentration aufrufen

- 3. Vergrößerten Ausschnitt anzeigen:
- Dazu wird im gewünschten Bereich des Diagramms mit dem Finger/Stift ein Rechteckbereich aufgezogen. Die Größe des Rechtecks wird über das Ziehen einer Diagonalen von Startpunkt (Druckpunkt, links am oberen Rahmen des Diagramms) und Endpunkt (Loslassen, rechts am unteren Rahmen des Diagramms) bestimmt.
- Auf eine beliebige Stelle innerhalb des markierten Rechteckbereichs drücken. Der Ausschnitt wird dann vergrößert angezeigt.
- Der Vorgang kann beliebig oft wiederholt werden, bis der Ausschnitt in der gewünschten Vergrößerung angezeigt wird oder die maximale Vergrößerungsstufe (max. 30 Datenlogger-Elemente, entspricht 30 min Verlauf bei einem Speicherzyklus von 60 s) erreicht ist.
- Im Zoom-Modus kann zusätzlich vor- und zurück geblättert werden.

- 4. Wieder den Gesamtverlauf anzeigen:
- Rechteck über einen kleinen Bereich des Diagramms aufziehen und auf eine beliebige Stelle außerhalb des markierten Bereichs drücken.
- 5. Verlaufsanzeige schließen:
- Taste **Ende** drücken.
- Es erfolgt ein Rücksprung ins Hauptmenü.

Hinweis Speicherzyklus der Aufzeichnung:

Das Zeitintervall der Datenaufzeichnung kann im Dialogfenster **Speicherzykluszeit** ("Speicherzyklus einstellen:" auf Seite 6-29) neu definiert werden.

Fehlermeldungen

Die Fehlerdetektion ist Teil des geräteinternen Kontrollsystems. Es überwacht die Regelkreise inklusive deren Sensorik. Wird ein Fehler im System ermittelt, wird das Alarmrelais geschaltet und löst folgende Signal- und Meldevorgänge aus:

- ein akustischer Alarm wird als abgesetzter Hupton ausgegeben,
- ein blinkendes Warndreieck und das entsprechende Icon werden am Hauptmenü eingeblendet; die Wertanzeigen werden dann nicht mehr aktualisiert,
- der detektierte Fehler wird in der Fehlertabelle gelistet,
- der Vorgang wird in die Ereignisanzeige eingetragen.
- Steht ein Fehler, nachdem er bestätigt wurde, weiterhin an, kann dieser Fehler über eine rote Taste (T, CO₂, O₂, RH oder System, Menütaste) wieder ausgelesen werden.

Reaktion auf das Ereignis Fehlermeldung

Wurde das Alarmrelais aufgrund einer Bedienaktion geschaltet, kann der Schaltzustand durch Bestätigung der Fehlermeldung zurückgesetzt werden (z.B. beim manuellen Abbruch der Steri-run-Dekontaminationsroutine).

Wurde das Alarmrelais aufgrund eines technischen Fehlers geschaltet, bleibt der Schaltzustand solange aktiviert, bis die Fehlerursache beseitigt ist (z.B. Wasserstand in Nutzraum zu niedrig).

- 1. Fehlermeldung bestätigen:
- Wird das Warndreieck eingeblendet, auf eine beliebige Stelle des Tastbildschirms drücken.
- Das Dialogfenster Fehler wird angezeigt und die detektierte Fehlerursache eingeblendet.
- Das akustische Signal wird abgeschaltet.
- 2. Fehleranzeige schließen:
- Taste Ende drücken.
- Fehlermeldung wird ausgeblendet.



Abbildung 6-51. Ereignis Fehlermeldung

Übertemperaturschutz zurücksetzen



Hat die Gerätesteuerung den Übertemperaturschutz aktiviert und auf Notregelung umgeschaltet, werden ein blinkendes Warndreieck und das Icon im Hauptmenü eingeblendet.



Abbildung 6-52. Fehlermeldung Übertemperatur

- 1. Fehlerursache anzeigen:
- Auf eine beliebige Stelle des Tastbildschirms drücken.
- Das Dialogfenster Fehler wird angezeigt und die detektierte Fehlerursache eingeblendet.
- Das akustische Signal wird abgeschaltet.

Fehler	
Temp - Istwert über	
-	
Ende	

Abbildung 6-53. Fehlermeldung Übertemperatur

- 2. Fehleranzeige schließen:
- Taste Ende drücken.
- Fehlermeldung wird ausgeblendet.
- Temperatur-Anzeigefeld wird rot umrandet.
- 3. Fehlermeldung zurücksetzen:
- Gerät ausschalten.
- 4. Türen öffnen und Nutzraum abkühlen lassen.
- 5. Gerät einschalten.

Wird der Übertemperaturschutz trotz der Beseitigung möglicher Fehlerursachen (siehe Fehlertabelle) erneut aktiviert, das Gerät außer Betrieb nehmen und den technischen Service anfordern.

Maßnahmen nach einem Stromausfall

Nach Ausfall der Spannungsversorgung kann es zur Kondensation der Innenraumfeuchte an den Sensoren kommen. Dies kann ihre Funktion so weit beeinträchtigen, dass falsche Messwerte angezeigt werden oder das Gerät sogar einen Defekt meldet (Fühlerbruch, siehe "Übersicht Fehlerursachen und Fehlerbeseitigung" auf Seite 6-48).

Um einen fehlerfreien Betrieb zu gewährleisten, müssen folgende Maßnahmen getroffen werden:

- 1. Wasser ablassen und Innenraum abtrocknen.
- 2. Gerät ohne Wasser für 1 Stunde auf 55 °C aufheizen lassen.
- 3. Danach Gerät mit geöffneten Türen abkühlen lassen.
- 4. Anschließend Inbetriebnahme auf Inkubationstemperatur gemäß Kapitel "Inbetriebnahme" auf Seite 4-1.

Alternativ dazu oder wenn die vorstehend beschriebenen Maßnahmen nicht den gewünschten Erfolg bringen, kann auch ein Desinfektionsvorgang bei 180 °C gestartet werden. Siehe hierzu Kapitel "Steri-run-Dekontaminationsroutine" auf Seite 8-7.

Der Desinfektionsvorgang kann nach ungefähr 1 Stunde abgebrochen werden. Die Sensoren sollten bis dahin wieder trocken sein.

Übersicht Fehlerursachen und Fehlerbeseitigung

Die Fehlertabellen geben Auskunft über Fehlerquelle, Fehlerursache und mögliche Fehlerbeseitigung. Zur Kommunikation mit dem Technischen Service von Thermo Fisher Scientific halten Sie bitte die Gerätedaten bereit.

Regel- kreis	Fehlermeldung	Ursache	Abhilfe	Alarm- relais	Hupe	Log
System	Gerätetür zu lange offen	Gerätetür ist länger als 10 Min. geöffnet	Gerätetür schließen	Х	Х	Х
	Fehler: Display	Display kommuniziert nicht mit dem Mainboard ^{*1)}	Gerätereset. Bei wiederholtem Auftreten den Service anfordern	Х	Х	Х
	Fehler: EEPROM Mainboard	EEPROM auf Mainboard ist defekt	Gerätereset. Bei wiederholtem Auftreten den Service anfordern	Х	Х	Х
	Fehler: Datenlogger	Fehler beim Schreiben auf den Speicher des Datenloggers. Inkubator ist weiter funktionsfähig.	Gerätereset. Bei wiederholtem Auftreten den Service anfordern			
	Fehler: steri-run	Fehler in der Steri-run-Routine	Gerätereset. Bei wiederholtem Auftreten den Service anfordern	Х	Х	Х
	Power-Down während steri-run	Stromausfall während der Steri-Run Routine	Gerät neu starten und Steri-run erneut ausführen.	Х	Х	Х
	Fehler: auto-start	Fehler in der auto-start-Routine	Auto-start erneut ausführen. Bei wiederholtem Auftreten den Service anfordern.	Х	Х	Х
	Fehler: ADC	Messung des Referenzwiderstandes außerhalb der Toleranz	Gerätereset. Bei wiederholtem Auftreten den Service anfordern.	Х	Х	Х
	Fehler: Lüfter	Istwert des Lüfters liegt außerhalb der Toleranz.	Gerätereset. Bei wiederholtem Auftreten den Service anfordern.	Х	Х	Х
	IR-Sensor gewechselt	Neue Seriennummer erkannt	Alarm quittieren	Х	Х	Х

Regel- kreis	Fehlermeldung	Ursache	Abhilfe	Alarm- relais	Hupe	Log
Tempera- tur	Sensor-/ Fühlerbruch	Messwert liegt außerhalb des akzeptierten Limits	Service anfordern. Trocknungsprozess durchführen.	Х	Х	Х
	Istwert über	Istwert > Sollwert + 1°C *2) *4)	Zulässige Umge- bungstemperatur nicht überschreiten / Service anfordern.	Х	Х	Х
	Istwert unter	Istwert < Sollwert + 1 °C ^{*3) *4)}	Setzt sich der Fehler nicht automatisch zurück, den Service anfordern.	Х	Х	Х
	Istwert nicht plausibel	Plausibilität des Temperatursignals nicht mehr gegeben	Gerätereset. Bei wie- derholtem Auftreten den Service anfordern.	Х	Х	Х
	Kalibrierwerte zu groß/klein	Max. Abgleichwert für Temperatur über- / unterschritten	Alarm quittieren, anderen Zielwert eingeben.			Х

Regel- kreis	Fehlermeldung	Ursache	Abhilfe	Alarm- relais	Hupe	Log
020% CO ₂	Sensor-/ Fühlerbruch	Messwert liegt außerhalb des akzeptierten Limits	Auto-start durchfüh- ren. Wenn im Anschluss der Fehler erneut auftritt, Feh- lerbehebung gemäß Kapitel "Maßnah- men nach einem Stromausfall" auf Seite 6-47 durchfüh- ren. Wenn sich der Fehler nicht beheben lässt, den Service anfordern.	Χ	Χ	Χ
	Istwert über	Istwert > Sollwert + 1% *4)	automatisch	Х	Х	Х
	Istwert unter	Istwert < Sollwert - 1% *3) *4)	automatisch	Х	Х	Х
	RH Störung Kommunikation	RH Sensor kommuniziert nicht mit dem Mainboard	automatisch	Х	Х	Х
	Kalibrierwerte zu groß/klein	Max. Abgleichswert für CO ₂ über- / unterschritten	Alarm quittieren			Х
	Störung Kommunikation	Sensor kommuniziert nicht mit dem Mainboard	automatisch	Х	Х	Х
	Fehler: Gasflaschenumschalter	Gasflaschenumschalter kommuniziert nicht mit dem Mainboard	automatisch	Х	Х	Х
	Kein Gas	Beide CO ₂ -Gasflaschen sind leer	Mindestens eine CO ₂ -Flasche tauschen.	Х	Х	Х
	RH Sensor-/Fühlerbruch	Messwert liegt außerhalb des akzeptierten Limits	Service anfordern. Siehe auch Kapitel "Maßnahmen nach einem Stromausfall" auf Seite 6-47.	Х	Х	Х
	Gasflasche A leer	Gasflasche A ist leer	Gasflasche A wechseln			
	Gasflasche B leer	Gasflasche B ist leer	Gasflasche B wechseln			

Regel- kreis	Fehlermeldung	Ursache	Abhilfe	Alarm- relais	Hupe	Log
0100% O ₂	Sensor-/ Fühlerbruch	Messwert liegt außerhalb des akzeptierten Limits	Service anfordern	Х	Х	Х
	Istwert über	Istwert > Sollwert + 1% *4)	Gasversorgung überprüfen. Vordruck auf max. 1 bar verringern.	Х	Х	Х
	Istwert unter	Istwert < Sollwert - 1% *4)	Gasversorgung überprüfen. Gasflasche tauschen. Vordruck auf max. 1 bar vergrößern. Zuleitung überprüfen.	Χ	Х	Х
	Störung Kommunikation	Sensor kommuniziert nicht mit dem Mainboard	Service anfordern	Х	Х	Х
	Fehler: Gasflaschenumschalter	Gasflaschenumschalter kommuniziert nicht mit dem Mainboard	automatisch	Х	Х	Х
	Kein Gas	Beide O ₂ -Gasflaschen sind leer	Mindestens eine O ₂ -Flasche tauschen.	Х	Х	Х
	Gasflasche A leer	Gasflasche A ist leer	Gasflasche A wechseln			
	Gasflasche B leer	Gasflasche B ist leer	Gasflasche B wechseln			

Regel- kreis	Fehlermeldung	Ursache	Abhilfe	Alarm- relais	Hupe	Log
rH	Kein Wasser	Zu wenig Wasser im Wasserreservoir.	Wasser nachfüllen oder, wenn trockener Betrieb gewünscht ist, den Wasserstandsensor deaktivieren. Wenn die Fehlermeldung trotz Nachfüllen wiederkehrt, Fehlerbehebung gemäß Kapitel "Maßnahmen nach einem Stromausfall" auf Seite 6-47 durchführen. Wenn sich der Fehler nicht beheben lässt, den Service anfordern.	x	X	X

^{*1)} Der Fehler wird nur über das Display ausgegeben und nicht in die Fehlertabelle eingetragen.

*2) Bei Auftreten des Fehlers wird eine Spezialregelung zum Schutz der Proben aktiv. Um dies zu verdeutlichen erscheint das Icon.

*3) Wartezeit bis zur Fehlermeldung:

- 45 Min. nach einer Türöffnung,

- 159 Min. nach einer Sollwertänderung.

*4) Dieser Wert kann durch den Service verändert werden.

Außerbetriebnahme

Inhalt

• "Gerät außer Betrieb nehmen" auf Seite 7-1

Gerät außer Betrieb nehmen



WARNUNG Kontaminationsgefahr! Die Oberflächen des Nutzraumes können kontaminiert sein. Es besteht die Gefahr, dass Keime auf die Umgebung übertragen werden. Gerät zur Außerbetriebnahme dekontaminieren!

- 1. Kulturgefäße mit den Kulturen und alle Hilfsmittel aus dem Nutzraum herausnehmen.
- 2. Auffanggefäß mit ausreichendem Fassungsvermögen bereitstellen.
- 3. Schlauchende in das Auffanggefäß halten und Schlauchventil sicher einrasten.
- 4. Das Wasserreservoir beginnt sich zu entleeren.



Abbildung 7-1. Befüll- und Ablassventil des Wasserreservoirs

- 5. Wasserreservoir vollständig in das Auffanggefäß entleeren lassen.
- 6. Schlauchventil wieder entfernen.
- 7. Steri-run-Dekontaminationsroutine starten ("Steri-run aufrufen" auf Seite 6-16).

- 8. Gerät am Netzschalter abschalten, wenn die Steri-run-Dekontaminationsroutine beendet ist.
- 9. Netzstecker aus der Steckdose ziehen und gegen Wiederanschluss sichern.
- 10. Schließventile der CO $_2$ /O $_2$ /N $_2$ -Versorgungsanlage schließen.
- 11. Gasdruckschläuche vom Anschlussventil an der Geräterückseite abziehen.
- 12. Während der Zeitphase der Stilllegung des Gerätes muss der Nutzraum ständig belüftet werden. Dazu die Glastür und die Außentür leicht öffnen und im geöffneten Zustand sichern.

8

Reinigung und Desinfektion

Inhalt

- "Reinigung" auf Seite 8-1
- "Dekontaminationsverfahren" auf Seite 8-2
- "Desinfektion oder Steri-run vorbereiten" auf Seite 8-2
- "Wisch-/Sprühdesinfektion" auf Seite 8-3
- "Steri-run-Dekontaminationsroutine" auf Seite 8-7

Reinigung

VORSICHT Unverträgliche Reinigungsmittel!



Teile des Gerätes sind aus Kunststoffen gefertigt. Lösemittel können Kunststoffe anlösen. Starke Säuren oder Laugen können eine Versprödung der Kunststoffe verursachen. Zum Reinigen der Kunststoffteile und -oberflächen keine kohlenwasserstoffhaltigen Lösemittel, keine Mittel mit einem Alkoholgehalt von mehr als 10% und keine starken Säuren oder Laugen benutzen!

Feuchteempfindliche Bauteile!

Das Display sowie den Schaltkasten an der Geräterückseite nicht mit Reinigungsmittel besprühen. Beim Abwischen darauf achten, dass keine Feuchtigkeit in diese Bauteile eindringt.

Reinigung der Außenflächen:

- 1. Schmutzrückstände und Ablagerungen mit lauwarmem Wasser, das mit handelsüblichem Spülmittel versetzt ist, gründlich beseitigen.
- 2. Die Oberflächen mit einem sauberen Tuch und klarem Wasser abwischen.
- 3. Abschließend die Oberflächen mit einem sauberen Tuch trocken reiben.

Reinigung des Displays:



VORSICHT Feuchteempfindliches Display! Das Display nicht feucht abwischen oder mit Reinigungsmittel besprühen!

• Das Display mit einem Tuch aus 100 % Microfaser trocken abwischen!

Dekontaminationsverfahren

Informationen zur Reinigung und Desinfektion des Cell Locker sind in der separaten Betriebsanleitung des Cell Locker zu finden (siehe Anhang).

Zur Dekontamination des Gerätes müssen vom Betreiber Hygienerichtlinien bekanntgegeben werden, die die Dekontaminationsmaßnahmen auf den Einsatz des Gerätes abstimmen. Folgende Desinfektionsverfahren sind für das Gerät geeignet:

Wisch-/Sprühdesinfektion:

Die Wisch-/Sprühdesinfektion ist für das Gerät und für alle Hilfsmittel als standardisiertes manuelles Desinfektionsverfahren vorgesehen.

Die Steri-run-Dekontaminationsroutine dekontaminiert in einem automatisierten Programmablauf den kompletten Nutzraum inklusive des Regalsystems und der Sensoren.

Desinfektion oder Steri-run vorbereiten

- 1. Sämtliche Proben aus dem Probenraum herausnehmen und sicher einlagern.
- 2. Auffanggefäß mit ausreichendem Fassungsvermögen bereitstellen.
- 3. Schlauchende in das Auffanggefäß legen und das Schlauchventil sicher einrasten.
- 4. Das Wasserreservoir beginnt sich zu entleeren.



Abbildung 8-1. Befüll- und Ablassventil des Wasserreservoirs

- 5. Wasserreservoir vollständig in das Auffanggefäß entleeren lassen.
- 6. Wasserrückstände mit einem Tuch aufwischen.
- 7. Airbox (1/Abbildung 8-2) vom Sockel der Wasserreservoirabdeckung abziehen und herausnehmen.
- 8. HEPA-Filter (2/Abbildung 8-2) aus Airbox (1/Abbildung 8-2) ausbauen und Airbox wieder einbauen.

Wisch-/Sprühdesinfektion

Die manuelle Wisch-/Sprühdesinfektion wird in drei Arbeitsabschnitten durchgeführt:

- Vordesinfektion
- Reinigung
- Enddesinfektion

VORSICHT

 Alkoholische Desinfektionsmittel! Desinfektionsmittel, die mehr als 10 % Alkohol enthalten, können zusammen mit Luft leicht entflammbare und explosive Gasgemische bilden.
Bei Anwendung solcher Desinfektionsmittel offenes Feuer oder starke Hitzeeinwirkung

Bei Anwendung solcher Desinfektionsmittel offenes Feuer oder starke Hitzeeinwirkung während des gesamten Desinfektionsverfahrens vermeiden!

- Desinfektionsmittel ausschließlich in gut belüfteten Räumen anwenden.
- Nach Einwirkung des Desinfektionsmittels die behandelten Geräteteile gut trocken reiben.
- Sicherheitsregeln zur Vermeidung von Brand- und Explosionsgefahren durch alkoholische Desinfektionsmittel beachten.



VORSICHT

Chloridhaltige Mittel! Keine chloridhaltigen Desinfektionsmittel verwenden!



WARNUNG Stromschlag!

Die Berührung stromführender Teile kann zu einem lebensgefährlichen Stromschlag führen. Vor Beginn der manuellen Reinigungs- und Desinfektionsarbeiten das Gerät vom Netz nehmen!

- Gerät am Netzschalter abschalten.
- Netzstecker ziehen und gegen versehentlichen Wiederanschluss sichern.
- Kontrollieren, ob das Gerät spannungsfrei ist.



VORSICHT Gesundheitsgefährdung!

Die Oberflächen des Nutzraums können kontaminiert sein. Der Kontakt mit kontaminierten Reinigungsflüssigkeiten kann Infektionen verursachen. Desinfektionsmittel können gesundheitsschädliche Stoffe enthalten.

Bei der Reinigung und Desinfektion die Schutzmaßnahmen und Hygieneregeln einhalten!

- Schutzhandschuhe tragen.
- Schutzbrille tragen.
- Zum Schutz der Schleimhäute Mund- und Nasenschutz tragen.
- Beachten Sie die Hinweise des Desinfektionsmittelherstellers und der Hygienefachkraft.

Vordesinfektion:

1. Die Oberflächen des Probenraumes und der Einbauten mit Desinfektionsmittel besprühen und abwischen.



VORSICHT Feuchtigkeitsempfindliche Bauteile!

Den CO₂-Sensor und den O₂/N₂Sensor hinter der Luftführung nicht mit Desinfektionsmittel besprühen.

2. Desinfektionsmittel entsprechend den Herstellerangaben einwirken lassen.

Einbauten und Regalsystem ausbauen:

- 1. Die Einlagebleche herausnehmen und anschließend das komplette Regalsystem aus dem Nutzraum ausbauen. Ein- und Ausbau des Regalsystems sind im Abschnitt "Regalsystem installieren" auf Seite 4-9 beschrieben.
- 2. Airbox und HEPA-Filter ausbauen. Ein- und Ausbau der Airbox und des HEPA-Filters sind im Abschnitt "HEPA-Filter tauschen" auf Seite 9-7 beschrieben.


Abbildung 8-2. HEPA-Filter und Airbox

- 3. Oberteil der Luftführung (1/Abbildung 8-3) in Richtung Gerätevorderseite ziehen und nach unten wegheben, wenn die Schlüssellochprofile an den vorderen Laschen die Haltestifte in der Nutzraumdecke freigeben.
- 4. Oberteil aus Rückteil der Luftführung (2/Abbildung 8-3) aushaken und aus dem Nutzraum herausnehmen.





Abbildung 8-3. Luftführung

- 5. Rückteil der Luftführung (2/Abbildung 8-3) aus der Rückwand aushängen und herausnehmen.
- 6. Vorfilter (9/Abbildung 3-9) herausnehmen. Der Vorfilter kann autoklaviert werden.
- 7. Abdeckung des Wasserreservoirs abbauen.

Reinigung des Nutzraumes und der Ausbauteile:

- 1. Schmutzrückstände und Ablagerungen mit lauwarmem Wasser, das mit handelsüblichem Spülmittel versetzt ist, gründlich beseitigen.
- 2. Die Oberflächen mit einem sauberen Tuch mit reichlich klarem Wasser abwischen.
- 3. Anschließend das Reinigungswasser aus dem Wasserreservoir entfernen und sämtliche Oberflächen des Nutzraumes gut trockenreiben.
- 4. Ausbauteile reinigen und ebenfalls gut trockenreiben.

Enddesinfektion:

- 1. Die Oberflächen des Nutzraumes, das Regalsystem und Ausbauteile erneut mit Desinfektionsmittel besprühen bzw. abwischen.
- 2. Desinfektionsmittel entsprechend den Herstellerangaben einwirken lassen.
- 3. Regalsystem und Ausbauteile wieder einbauen.

Steri-run-Dekontaminationsroutine

Steri-run ist eine automatisch ablaufende Sterilisationsroutine mit fester Nenntemperatur und einem softwaregesteuerten Ablauf für das Aufheizen, Halten auf Nenntemperatur und Abkühlen.

Der gesamte Programmdurchlauf der Dekontaminationsroutine dauert weniger als 12 Stunden. Während dieser Routine wird im Nutzraum für 90 Minuten eine trocken-heiße Atmosphäre bei 180 °C mit hoher dekontaminierender Wirkung erzeugt. Die Wirksamkeit der Steri-run-Dekontaminationsroutine wurde von unabhängigen Instituten nachgewiesen. Es wird eine Keimreduzierung um 10⁶ (6-log-Reduzierung) gemäß Norm ISO 11138 erreicht. Thermo Scientific stellt Informationen zu diesen Tests auf Anfrage zur Verfügung.

Der elektromechanische Türverschluss bietet bei der steri-run-Dekontaminationsroutine einen zusätzlichen Schutz gegen Verbrennungen. Er verriegelt die Außentür, wenn die Nutzraumtemperatur 65 °C erreicht, hält diese verschlossen, solange im Nutzraum gefährliche Temperaturen vorherrschen und entriegelt die Tür schließlich wieder, wenn die Nutzraumtemperatur wieder unter 65 °C sinkt.

Nach abgeschlossenem Durchlauf muss das Gerät mit der auto-start-Routine wieder in Betrieb genommen werden.

Hinweis Verhinderung des Starts der Steri-run-Dekontaminationsroutine: Die Steri-run-Dekontaminationsroutine kann nicht gestartet werden, wenn einer der folgenden Fehler detektiert ist.

Regelkreis Temperatur:

- Sensor-/Fühlerbruch,
- Außen- und/oder Innentür (bei offener Außentür erscheint die Fehlermeldung "Tür geöffnet"),
- Istwert über (Abweichung vom Sollwert zu stark),
- Istwert unter (Abweichung vom Sollwert zu stark),
- Istwert nicht plausibel,
- Störung Kommunikation,
- Wasser wurde detektiert.

Übertemperaturschutz:

Wurde bei dem Gerät der Übertemperaturschutz aktiviert, kann die Steri-run-Dekontaminationsroutine erst gestartet werden, nachdem der Fehler behoben bzw. zurückgesetzt wurde.

Keine Gasversorgung während der Steri-run-Dekontaminationsroutine bei Ausstattung mit Gasmonitoring-System (optional):

Tritt während der Steri-run-Dekontaminationsroutine der Fehler "Kein Gas" auf, wird das akustische Signal (Hupe) aktiviert. Das Signal kann durch Drücken auf eine beliebige Stelle des Displays bestätigt werden. In diesem Fall wird die Steri-run-Dekontaminationsroutine nicht abgebrochen. Das Alarmrelais bleibt so lange geschaltet, bis der Gasmonitor wieder eine volle Flasche erkennt.



VORSICHT Betriebstemperatur-Grenzwerte des Cell Lockers! Vor der Durchführung eines Dekontaminationslaufs sind die Cell Locker zu entfernen. Die max. Betriebstemperatur beträgt 121°C/250 °F.

Ablauf einer Steri-Run-Dekontaminationsroutine:

- 1. Vor dem Start der Dekontaminationsroutine entfernen Sie den Silikonstopfen vom Innenbehälter und stecken ihn in die Außenseite der Zugangsöffnung.
- 2. Die Komponenten des Regalsystems nach der Reinigung wieder in den Nutzraum einbauen.
- 3. Gerät am Netzschalter einschalten.
- 4. Dekontaminationsroutine aktivieren und starten.
- 5. Nach Ablauf der Steri-run-Routine das Gerät abschalten.
- 6. Airbox (1/Abbildung 8-2) ausbauen und HEPA-Filter wieder einbauen (2/Abbildung 8-2).
- 7. Bei Bedarf mit auto-start den Betrieb wieder aufnehmen.

VORSICHT Heiße Oberflächen!



Der Griff und die Scheibe der Glastür, das Innenblech der Außentür sowie die Oberflächen des Regalsystems und des Nutzraumes, werden während der Steri-run-Dekontaminationsroutine stark erhitzt.

Während des Ablaufes oder unmittelbar nach Abbruch der Routine diese Oberflächen nur mit Schutzhandschuhen anfassen!

Inkubatoren mit Türverschluss-Kit verhindern den Zugriff auf den Nutzraum, solange die Oberflächen im Innenraum durch die Steri-run-Dekontaminationsroutine heiß sind.

VORSICHT Beschädigung der Proben!



Der Nutzraum wird während der Steri-run-Dekontaminationsroutine auf 180 °C aufgeheizt. Sicherstellen, dass:

- alle Proben aus dem Nutzraum entnommen wurden,
- alle Hilfsmittel aus dem Nutzraum entfernt wurden.

Betriebsphasen der Steri-run-Dekontaminationsroutine:

Die Restlaufzeit der Steri-run-Dekontaminationsroutine bezeichnet die Zeitspanne vom Start bzw. dem aktuellen Zeitstatus der Routine bis zum Ende der Phase Abkühlen. Die angezeigten Restlaufzeiten sind keine Messwerte, sondern lediglich Orientierungswerte. Die Routine ist in drei Phasen gegliedert:

- 1. Anheizphase,
- 2. Dekontaminationsphase,
- 3. Abkühlen.

Anheizphase: ca. 2 Std.

Der Nutzraum wird auf 180 °C aufgeheizt. Inkubatoren, die mit dem elektromagnetischen Türverschluss-Kit ausgestattet sind, verriegeln die Tür, wenn die Nutzraumtemperatur 65 °C übersteigt.

Dekontaminationsphase: ca. 1.5 Std. Nach Erzeugung der Dekontaminationstemperatur wird die Dekontaminationsphase für ca. 90 Minuten gestartet. Dabei wird die Temperatur von 180 °C gehalten. Abkühlphase: ca. 8 Std.

Das Gerät kühlt wieder auf den ursprünglich eingestellten Temperatur-Sollwert ab. Inkubatoren, die mit dem elektromagnetischen Türverschluss-Kit ausgestattet sind, entriegeln die Türverriegelung, wenn die Nutzraumtemperatur unter 65°C sinkt.



Abbildung 8-4. Phasen der Dekontaminationsroutine

Steri-run aktivieren

Steri-run ist eine automatische Dekontaminationsroutine zur Desinfektion des Gerätenutzraumes.

1. Taste **steri-run** drücken.

Hinweis

Um Schäden am Silikon zu vermeiden, entfernen Sie unbedingt den Silikonstopfen aus dem Innenbehälter und stecken Sie ihn in die Außenseite der Zugangsöffnung, bevor Sie mit der Steri-run-Dekontamination beginnen.

• Das Menü steri-Run-Ablauf wird angezeigt.



Abbildung 8-5. Menü steri-run - Ablauf

- 2. Menü steri-run Ablauf verlassen und steri-run abbrechen:
- Taste **Ende** drücken.
- Es erfolgt ein Rücksprung ins Hauptmenü.

- 3. Steri-run aktivieren:
- Die Taste START betätigen.
- Das Dialogfenster Steri-Run-Ablauf wird angezeigt.
- 4. Nutzraum belüften, beide Gerätetüren öffnen, bis nach 30 s das Zeitsignal ertönt.
- 5. Alle Proben aus dem Nutzraum nehmen.
- 6. Das Wasser aus dem Wasserreservoir ablassen und Wasserrückstände aufwischen.
- 7. Nach Ertönen des Zeitsignals beide Gerätetüren schließen.
- Steri-Run starten.
- Der Ablauf der Steri-Run-Dekontaminationsroutine beginnt. Das Gerät heizt sich auf und der elektromechanische Türverschluss wird bei 65 °C aktiviert.
- Während des Verlaufs der Steri-run-Dekontaminationsroutine wird am Display der aktuelle Status angezeigt und dazu folgende Informationen ausgegeben:
 - Temperatur,
 - Startzeit,
 - Phase,
 - Restlaufzeit.



Abbildung 8-6. Steri-run aktivieren

Steri-run abbrechen

Die Steri-run-Dekontaminationsroutine kann jederzeit abgebrochen werden.

- 1. Steri-run abbrechen:
- Taste Stop drücken.
 - Wurde die Taste **Stop** gedrückt, wird als Sicherheitsabfrage das Dialogfenster steri-run Stop aufgerufen. Die Routine kann jetzt endgültig abgebrochen oder wieder fortgesetzt werden.
- 2. Steri-run beenden:
- Taste Ende drücken.
- Die Fehlermeldung wird angezeigt.
- Wird die Fehlermeldung bestätigt, erfolgt ein Rücksprung ins Hauptmenü.

- 3. Steri-run fortsetzen:
- Taste **Zurück** drücken.
- Es erfolgt ein Rücksprung in die Statusanzeige, die Dekontaminationsroutine wird fortgesetzt.
- 4. Steri-run aus der Statusanzeige heraus unterbrechen:
- Taste **Stop** drücken.
- Als Sicherheitsabfrage wird das Dialogfenster steri-run Stop aufgerufen. Weiter mit Arbeitsschritt 2.

Steri-run Fehlerabbruch

Tritt während es Ablaufs der Dekontaminationsroutine ein Fehler auf, wird eine Fehlermeldung ausgegeben und folgende Aktionen eingeleitet:

- Die Dekontaminationsroutine wechselt automatisch in die Kühlphase.
- Das akustische Signal (Hupe) wird ausgegeben.
- 1. Akustisches Signal bestätigen:
- Display an einer beliebigen Stelle drücken.
- Das akustische Signal verstummt. Die Taste **Ende** wird eingeblendet. Wird im Anschluss die Dekontaminationsroutine nicht manuell abgebrochen, wird auf die Solltemperatur abgekühlt und der Nutzraum getrocknet.
- 2. Steri-run abbrechen:
- Taste **Ende** drücken.
- Die Fehlermeldung wird angezeigt.
- Wird die Fehlermeldung bestätigt, erfolgt ein Rücksprung ins Hauptmenü.

steri-ru	n - At	bruch		
<ende></ende>	Stoppe	n der Desinfe	ktions-	
<zurück></zurück>	Fortset	zen		
		0		
	_		e	

Abbildung 8-7. Steri-run unterbrechen, abbrechen

Steri-run abschließen

Nach vollständigem Ablauf der 3 Phasen wird automatisch das Dialogfenster **steri-run - Ende** (Abbildung 8-8) angezeigt. Die Dekontaminationsroutine muss manuell beendet werden.

	steri-run - Ende
	1. Tür öffnen
	2. Wasser einfüllen
	3. Ggf. Einbauten wie HEPA-Filter einbauen 4. Tür schließen
	5. Ggf. auto-start-Routine durchführen
_	
	((Ende))

Abbildung 8-8. Steri-run beenden

- Steri-run beenden:
 - Taste Ende drücken.
 - Es erfolgt ein Rücksprung ins Hauptmenü.

Hinweis Tür öffnen während der steri-run-Dekontaminationsroutine:

Wird während der Dekontaminationsroutine die Tür geöffnet und anschließend wieder geschlossen, wird in diejenige Ablaufphase zurückgesprungen, die eine fehlerfreie Fortsetzung der Routine gewährleistet.

Hinweis

Verbrennungsgefahr! Das Öffnen der Tür ist, außer im Notfall, zu vermeiden.

Öffnen der Tür nach Abbruch der Steri-run-Dekontaminationsroutine

Bei Geräten mit elektromechanischem Türverschluss-Kit kann die Tür direkt nach dem Abbruch der steri-run-Dekontaminationsroutine durch den Benutzer oder aufgrund eines Fehlers nicht geöffnet werden.

Um die Verriegelung der Außentür zu entriegeln, bevor die Temperatur auf einen unbedenklichen Wert von 65 °C gesunken ist, ist der Notentriegelungshebel zu betätigen (Pos. 3 in Abbildung 8-9):

• Position des Notentriegelungshebels (3) an der Unterseite des Inkubators bestimmen. Er befindet sich an der linken Geräteseite, ca. 5 cm neben/hinter dem Netzschalter (4), (siehe Abbildung 8-9 unten).



Abbildung 8-9. Türverriegelungshebel und Notentriegelung an der Unterseite des Inkubators

- Unter die Bodenplatte greifen und den Notentriegelungshebels (3) ziehen.
- Der Türverschluss entriegelt und die Außentür kann geöffnet werden.

VORSICHT Heiße Oberflächen!



Der Griff und die Scheibe der Glastür, das Innenblech der Außentür sowie die Oberflächen des Regalsystems und des Nutzraumes, werden während der Steri-run-Dekontaminationsroutine stark erhitzt.

Während des Ablaufes oder unmittelbar nach Abbruch der Routine diese Oberflächen nur mit Schutzhandschuhen anfassen!

8 Reinigung und Desinfektion Steri-run-Dekontaminationsroutine

Instandhaltung

Inhalt

- "Inspektionen und Kontrollen" auf Seite 9-1
- "Wartungsintervalle" auf Seite 9-2
- "Temperaturabgleich vorbereiten" auf Seite 9-2
- "Temperaturabgleich durchführen" auf Seite 9-3
- "CO2-Abgleich vorbereiten" auf Seite 9-5
- "CO2-Abgleich durchführen" auf Seite 9-6
- "HEPA-Filter tauschen" auf Seite 9-7
- "Gaseinlassfilter tauschen" auf Seite 9-9
- "Austausch der Gerätesicherung" auf Seite 9-10
- "Türdichtung wechseln" auf Seite 9-10

Inspektionen und Kontrollen

Zur Erhaltung der Funktionstüchtigkeit und Sicherheit des Gerätes muss eine Prüfung der unten aufgeführten Funktionen und Gerätekomponenten in unterschiedlichen Zeitintervallen durchgeführt werden.

Informationen zur Instandhaltung des Cell Locker sind in der separaten Betriebsanleitung des Cell Locker zu finden (siehe Anhang).

Tägliche Kontrolle:

- Gasvorrat der CO₂-Versorgungsanlage.
- Gasvorrat der O₂-/N₂-Versorgungsanlage.

Jährliche Inspektion:

- Dichtigkeit der Glastürdichtung
- Durchlässigkeit der Druckausgleichsöffnung mit Einsatz
- Funktionstest des Bedienfeldes und der Geräteregelung
- Elektrische Sicherheitsprüfung entsprechend den national gültigen Vorschriften (z. B. BGV 3)

Hinweis Funktionsprüfung:

Wurden für Inspektionen Schutzeinrichtungen ausgebaut oder außer Funktion gesetzt, darf das Gerät erst wieder in Betrieb genommen werden, wenn die Schutzeinrichtungen wieder eingebaut und auf ihre einwandfreie Funktion hin überprüft wurden.

Wartungsintervalle

Im laufenden Betrieb sind folgende Wartungsarbeiten durchzuführen:

Vierteljährliche Wartung:

- auto-start-Routine und Steri-Run-Dekontaminationsroutine fahren.
- Temperatur- und CO₂-/O₂-Vergleichsmessung durchführen.

6-fach-Gasblende:

- Die Dichtung der Abdeckung alle 6 Monate prüfen und ggf. austauschen.
- Informationen zu den Membranfiltern des Cell Lockers sind in der entsprechenden Betriebsanleitung zu finden (siehe Anhang).

Jährliche Wartung:

- Gaseinlassfilter wechseln.
- Serviceprüfung des Technischen Services durchführen

Hinweis Wartungsvertrag:

Thermo Scientific bietet einen auf das Gerät abgestimmten Wartungsvertrag an, der alle erforderlichen Prüfungs- und Instandhaltungsdienste beinhaltet.

Temperaturabgleich vorbereiten

Zur Ermittlung des genauen Messwertes des geräteinternen Temperatursensors sollte vierteljährlich eine Temperaturvergleichsmessung durchgeführt werden. Wird dabei eine größere Messabweichung festgestellt, sollte ein Temperaturabgleich durchgeführt werden. Hierbei wird die Temperaturregelung des Gerätes auf den Messwert der Vergleichsmessung eingestellt.

Zur Vergleichsmessung sollte ein kalibriertes Messinstrument mit einer Genauigkeit < ± 0,1 °C verwendet werden.

Um zeitliche Temperaturschwankungen während der Messung zu minimieren, wird das Messinstrument in einem isothermen Behältnis (z. B. ein mit Glyzerin gefüllter Becher) im Nutzraum aufgestellt. Bezugsort der Vergleichsmessung ist die Mitte des Nutzraumes.

Hinweis Isothermes Behältnis:

Als isothermes Behältnis darf auf keinen Fall ein mit Wasser gefüllter Behälter verwendet werden, weil durch die Verdunstung von Wasser eine zu geringe Temperatur ermittelt wird.

Zu hohe Nutzraumtemperatur:

Eine möglicherweise zu hohe Nutzraumtemperatur nach dem Abgleich kann durch Öffnen der Türen für ca. 30 s abgebaut werden.



Abbildung 9-1. Temperaturabgleich vorbereiten

Vergleichsmessung durchführen:

- 1. Gerät am Netzschalter einschalten.
- 2. Temperatur-Sollwert einstellen und abwarten, bis das Gerät durchgewärmt ist. Dies kann mehrere Stunden dauern.
- 3. Messinstrument (2) mittig auf dem Einlageblech im mittleren Bereich des Nutzraumes aufstellen. Alternativ kann an gleicher Stelle ein Temperatursensor positioniert werden. Die Zuleitung wird entweder durch die Messöffnung in der Glastür (3) oder durch die Rohrdurchführung (1) an der Geräterückseite verlegt.
- 4. Türen schließen.
- 5. Abwarten, bis am Messinstrument ein konstanter Temperaturwert ablesbar ist.
- 6. Temperaturabgleich durchführen

Temperaturabgleich durchführen

Messbeispiel:

- Temperatur-Sollwert: 37 °C Gemessene Vergleichstemperatur: 36,4 °C
- 1. Taste Temperatur-Anzeigefeld drücken.
- Das Temperatur-Menü (Abbildung 9-2) wird angezeigt.



Abbildung 9-2. Temperatur-Anzeigefeld und Temperatur-Auswahlmenü

- 2. Temperatur-Menü wieder verlassen:
- Taste Ende drücken.
- 3. Untermenü Kalibrierung aufrufen:
- Taste Kalibration drücken.

Kalibration	Temperatur	kalibrierung	
	aktuell	37.0	
	neu	36.4	°C
	· ·	Enter	+

Abbildung 9-3. Temperaturabgleich durchführen

4. Gemessenen Wert (Zielwert) eingeben:

Der Zielwert kann stufenweise erhöht oder vermindert werden, durch Dauerdruck auf die jeweilige Taste - oder Taste + schaltet die Funktion in einen Schnelldurchlauf, nach ca. 3 s wird die Geschwindigkeit des Schnelldurchlaufes zusätzlich erhöht.

Zielwert erhöhen:

• Taste + drücken.

Sollwert verringern:

- Taste drücken.
- 5. Zielwert übernehmen und speichern:
- Taste Enter drücken.
- Taste **Speichern** drücken.

• Es erfolgt ein Rücksprung ins Hauptmenü. Der aktuelle im Nutzraum gemessene Istwert wird am Temperatur-Anzeigefeld angezeigt.

Hinweis Zu hohe Nutzraumtemperatur: Eine möglicherweise zu hohe Nutzraumtemperatur nach dem Abgleich kann durch Öffnen der Türen für ca. 30 s abgebaut werden.

Rücksetzung Wert:

Erfolgt 30 s lang keine weitere Änderung des Wertes, wird das Menü automatisch verlassen und der zuletzt bestätigte Wert wird übernommen.

CO₂-Abgleich vorbereiten

Zur Ermittlung eines genauen Messwertes des geräteinternen CO_2 -Sensors sollte vierteljährlich eine CO_2 -Vergleichsmessung durchgeführt werden.

Wird dabei eine größere Messabweichung festgestellt, sollte ein CO_2 -Abgleich durchgeführt werden. Hierbei wird die CO_2 -Regelung des Gerätes auf den Messwert der Vergleichsmessung eingestellt. Zur Vergleichsmessung sollte ein kalibriertes Messinstrument mit einer Genauigkeit von < \pm 0,3 % CO_2 verwendet werden.

Geeignetes Messgerät:

• Tragbares IR-Handmessgerät. Die Messprobe wird durch die verschließbare Messöffnung [1] der Glastür entnommen. Die Vergleichsmessung muss am durchgewärmten Gerät ausgeführt werden.

Vergleichsmessung durchführen:

- 1. Gerät am Netzschalter einschalten.
- 2. CO₂-Sollwert einstellen und warten, bis das Gerät durchgewärmt ist und sich die Feuchte aufgebaut hat.
- 3. Messsonde des IR-Handmessgerätes durch die Messöffnung in den Nutzraum führen. Abwarten, bis am Messgerät ein konstanter CO₂-Wert ablesbar ist.
- 4. Bei Geräten, die mit der optionalen Gasblende ausgerüstet sind, befindet sich die Messöffnung:
- bei STERI-CYCLE i160 in der mittleren Gasblende,
- bei **STERI-CYCLE i250** in der linken oder mittleren Gasblende.
- 5. Messsonde abziehen, die Messöffnung verschließen und die Türen schließen.
- 6. CO₂-Regelung kalibrieren.

Hinweis IR-Messzelle:

Bei Geräten mit IR-Messzelle kann der CO₂-Abgleich nur bei einer eingestellten CO₂-Konzentration von 4,0 % CO₂ oder höher durchgeführt werden. Die Kalibrierung sollte bei dem für den Arbeitsprozess vorgesehenen CO₂-Sollwert (künftiger Arbeitsprozesswert) vorgenommen werden.



Abbildung 9-4. Messöffnung Gasblenden

CO₂-Abgleich durchführen

Messbeispiel:

- CO₂-Sollwert: 5 % Referenzwert: 5,6 %
- 1. Taste CO₂-Anzeigefeld drücken.
- Das CO₂-Menü wird angezeigt.



- 2. CO₂-Menü wieder verlassen:
- Taste **Ende** drücken.
- 3. Untermenü Kalibrierung aufrufen:
- Taste Kalibrierung drücken.



Abbildung 9-5. CO 2-Abgleich durchführen

4. Gemessenen Wert (Zielwert) eingeben:

Der Zielwert kann stufenweise erhöht oder vermindert werden, durch Dauerdruck auf die jeweilige Taste - oder Taste + schaltet die Funktion in einen Schnelldurchlauf, nach ca. 3 s wird die Geschwindigkeit des Schnelldurchlaufes zusätzlich erhöht.

Zielwert erhöhen:

• Taste + drücken.

Sollwert verringern:

- Taste drücken.
- 5. Zielwert übernehmen und speichern:
- Taste Enter drücken.
- Taste Speichern drücken.
- Es erfolgt ein Rücksprung ins Hauptmenü. Der aktuelle, im Nutzraum gemessene Istwert wird am CO₂-Anzeigefeld angezeigt.

Hinweis Zu hoher CO₂-Gehalt:

Ein möglicherweise zu hoher CO₂-Gehalt nach dem Abgleich kann durch Öffnen der Türen für ca. 30 s abgebaut werden.

Rücksetzung Wert:

Erfolgt 30 s lang keine weitere Änderung des Wertes, wird das Menü automatisch verlassen und der zuletzt bestätigte Wert wird übernommen.

HEPA-Filter tauschen

Das HEPA-Filter sitzt unter einer Kunststoffhaube (Airbox) am Boden vor der Rückwand des Nutzraums.

Arbeitsschritte zum Austauschen des HEPA-Filters:

- 1. Gerät ausschalten, Gaszufuhr abstellen und Nutzraum ablüften lassen.
- 2. Mittleres und unteres Einlageblech aus Nutzraum herausnehmen.



3. Die Wasserreservoirabdeckung an der Frontseite anheben (1/Abbildung 9-6).

Abbildung 9-6. Airbox ausbauen

- 4. Die Wasserreservoirabdeckung nach vorn herausziehen (2/Abbildung 9-6).
- 5. Die Airbox mit HEPA-Filter abnehmen.
- 6. Airbox umdrehen und Laschen (5/Abbildung 9-7) auf der linken Seite der Airbox aus den Rasten (6/Abbildung 9-7) des HEPA-Filters aushaken.
- 7. Die Laschen auf der rechten Seite (3/Abbildung 9-7) der Airbox (1/Abbildung 9-7) aus den entsprechenden Nuten im HEPA-Filter (4/Abbildung 9-7) herausziehen.



Abbildung 9-7. HEPA-Filter einbauen

- 8. Neues HEPA-Filter in Airbox (1/Abbildung 9-7) einsetzen und einrasten lassen.
- 9. Airbox auf den Sockel der Wasserreservoirabdeckung aufsetzen.
- 10. Wurde das Gerät zuvor ohne HEPA-Filter betrieben, HEPA-Filter gemäß Anleitung "HEPA-Filter aktivieren / deaktivieren:" auf Seite 6-39 in der Benutzerkonfiguration aktivieren.
- 11. Ggf. das Erinnerungsintervall für den Austausch des HEPA-Filters gemäß Anleitung "Erinnerungsintervalle einstellen" auf Seite 6-26 in der Benutzerkonfiguration einstellen. Das Erinnerungsintervall für den Austausch des HEPA-Filters kann zwischen 1 und 12 Monaten eingestellt werden. Als Einstellwert wird die werkseitige Voreinstellung von 6 Monaten empfohlen.

Gaseinlassfilter tauschen

Das Gaseinlassfilter (CO_2 - $/O_2$ - $/N_2$ -Zufuhr) ist mit einem Kunststoffgewinde versehen und wird handfest in die Gewindeaufnahme am Schaltkasten geschraubt.

Arbeitsschritte für Gaseinlassfilter der Gaszufuhr:

- 1. Sicherstellen, dass die Gaszufuhr geschlossen ist.
- 2. Schlauchschelle (4/Abbildung 9-8) lösen.

3. Gasschlauch (5/Abbildung 9-8) von Anschlussstutzen des Gaseinlassfilters abziehen.

Arbeitsschritte für sämtliche Gaseinlassfilter:

- 4. Sicherungsblech (1/Abbildung 9-8) abschrauben.
- 5. Gaseinlassfilter (2/Abbildung 9-8 aus der Gewindeaufnahme (3/Abbildung 9-8) herausschrauben.
- 6. Beim Einschrauben des neuen Gaseinlassfilters darauf achten, dass das Kunststoffgewinde nicht verkantet angesetzt wird. Gaseinlassfilter vorsichtig einschrauben.
- 7. Sicherungsblech anschrauben.

Arbeitsschritte für Gaseinlassfilter der Gaszufuhr:

8. Gasschlauch auf den Anschlussstutzen des Filters setzen und mit der Schlauchschelle fixieren. Überprüfen, ob der Gasschlauch dicht auf dem Anschlussstutzen sitzt.



Abbildung 9-8. Gaseinlassfilter montieren

Austausch der Gerätesicherung

Ein Austausch der Gerätesicherungen durch den Anwender ist nicht möglich. Bei Ausfall des Gerätes infolge eines elektrischen Fehlers ist der Technische Service zu rufen.

Türdichtung wechseln

Hinweis

Es wird empfohlen, den Austausch der Türdichtung von einem Servicetechniker oder qualifiziertem Fachpersonal durchführen zu lassen.

Entsorgung

Inhalt

• "Übersicht der verwendeten Materialien:" auf Seite 10-1

WARNUNG Kontaminationsgefahr!



Das Gerät könnte zur Be- und Verarbeitung von infektiösen Substanzen eingesetzt worden sein. Das Gerät oder Teile des Gerätes können deshalb kontaminiert sein. Alle Komponenten des Gerätes müssen vor der Entsorgung dekontaminiert werden!

- Die Bauteile des Gerätes sind gründlich zu reinigen und anschließend abhängig vom Einsatzzweck entweder zu desinfizieren oder zu dekontaminieren.
- Dem Entsorgungsgut ist eine Unbedenklichkeitserklärung mit genauen Hinweisen über die Durchführung der Dekontaminationsmaßnahmen beizufügen.

Alle Komponenten des Gerätes können nach entsprechender Dekontaminierung der geregelten Entsorgung zugeführt werden.

Die HEPA-Filter müssen entsprechend den national gültigen Richtlinien für Sondermüll entsorgt werden.

Hinweis Recycling-Service: Thermo Fisher Scientific bietet einen umweltgerechten Recycling-Service für Altgeräte an.

Übersicht der verwendeten Materialien:

Komponente	Material
Thermische Isolationsteile	Glaswolle, Glaswolle mit einseitig kaschierten Glasvlies
Elektronikplatinen	Umhüllte elektrische Bauteile mit diversen Kunststoffen behaftet, auf epoxidharzgebundenen Leiterplatten bestückt.
Kunststoffteile, generell	ABS und PPS GF40, Materialkennzeichnung beachten

Komponente	Material
Außengehäuse	Stahlblech verzinkt und lackiert, Edelstahl 1.4016
Geräterückwand	Stahlblech verzinkt
Außentür	Stahlblech verzinkt und lackiert, Edelstahl 1.4016
Türinnenblech	Edelstahlblech 1.4301
Bedien- und Anzeigefolie	Polyethylen
Heizungen	Silikonummantelte Widerstandsheizleiter
Innenbehälter, Einbauten und Einlagebleche	Edelstahl 1.4301, Kupfer
Einsatz für Druckausgleichsöffnung	Edelstahl 1.4301 (Sockel), 1.4404 (Sinterfilter)
Glasscheibe	Natriumsilikatglas
Sensorblock (WLD)	Edelstahl 1.4301
Leitungen	Kunststoff- und Silikonummantelte Kupferlitze
Elastomere, generell	Silikon
Filter	HEPA-Filter, Glas (mikrofein), Cell Locker: Membranfilter mit ABS und Silikon Gasfilter Polypropylen Gehäuse und GF/PTFE-Membran, Vorfilter Metalldrahtgewebe Edelstahl 1.4401
Verpackung	Wellpappe, Polyethylenfolie, Styroporformteile und Polypropylen
Magnet der Türdichtung	Dauermagnet
Cell Locker	Polycarbonat Makrolon 2528

Technische Daten

Inhalt

- "STERI-CYCLE i160" auf Seite 11-2
- "STERI-CYCLE i250" auf Seite 11-6

STERI-CYCLE i160

Bezeichnung	Einheit	Wert
Mechanisch		
Außenabmessungen (B x H x T)	mm	637 x 905 x 790
Innenabmessungen (B x H x T)	mm	470 x 607 x 576
Innenraumvolumen, davon Nutzvolumen	1 1	ca. 165 ca. 100
Einlagebleche (B x T) Anzahl, Lieferumfang Anzahl, maximal Flächenlast, maximal	mm St. St. kg	423 x 465 3 11 10 / Einlageblech
Gesamtlast Gerät, maximal	kg	30
Gewicht	kg	83
Gewicht, ohne Zubehör	kg	78,4
Thermisch		
Thermische Sicherheitseinrichtungen nach DIN 12880:2007-05		Klasse 3.1 (Temperatur – Wählwächter (TWW) mit Wächterfunktion bei Temperaturüberschreitung)
Umgebungstemperaturbereich	°C	+1834
Umgebungstemperatur gestapelter Geräte	°C	+1828
Temperaturregelbereich, Inkubation	°C	RT + 355
Temperaturabweichung, zeitlich (DIN 12880, Teil 2) bei 37 °C	°C	± 0,1
Temperaturabweichung, räumlich (DIN 12880, Teil 2) bei 37 °C *1)	°C	<= ± 0,3
Dauer der auto-start-Routine: bis 37 °C Umgebungstemperatur 20 °C	Std.	510
Wärmeabgabe an Umgebung: bei 37 °C während Steri-Run-Dekontamination	kWh/h kWh/h	0,06 0,59
Feuchte		
Wasserqualitäten		Elektrischer Widerstand: 50 kOhmcm bis 1 MOhmcm Leitfähigkeit: 1 bis 20 μS/cm
Füllmengen: Inkubationsbetrieb	1	max. 3 / min. 0,5

Bezeichnung	Einheit	Wert
Konstantfeuchte bei 37 °C (High-Feuchte Modus) Konstantfeuchte bei 37 °C (Low-Feuchte Modus)	% rF % rF	ca. 93 ca. 90
Sonstiges		
Schalldruckpegel (DIN 45 635, Teil 1)	dB(A)	< 50
Relative Feuchte der Umgebung	% rF	max. 80
Höhe des Aufstellungsortes	m NN	max. 2000

^{*1)} Werte ermittelt in Anlehnung an DIN 12880 für Geräte in Standardausführung. Nähere Angaben sind der Kalibrieranweisung zu entnehmen

STERI-CYCLE i160

Bezeichnung	Einheit	Wert
CO ₂ -Gastechnik		
Gasreinheit	%	min. 99,5 oder med. Qualität
Vordruck	bar	min. 0,8 - max. 1
Mess- und Regelbereich	Vol -%	020
Regelabweichung, zeitlich	Vol -%	± 0,1
CO ₂ -Messzelle		
Genauigkeit bei 37 °C und 5% CO ₂	%CO ₂	± 0,3
O ₂ -Gastechnik		
Gasreinheit	%	min. 99,5 oder med. Qualität
Vordruck	bar	min. 0,8 - max. 1
Mess- und Regelbereich	Vol -%	1 21 bzw. 590
Regelabweichung, zeitlich	Vol -%	± 0,2
O ₂ -Messzelle		
Genauigkeit bei 37 °C und 21% O ₂	%O ₂	± 0,5 (Option: 121% O ₂) ± 2,0 (Option: 590% O ₂)
Elektrisch		
Nennspannung	V	1/N/PE 230 V, AC (± 10%) 1/N/PE 220 V, AC (± 10%) 1/N/PE 120 V, AC (± 10%) 1/N/PE 100 V, AC (± 10%)
Nennfrequenz	Hz	50/60
Schutzart (IEC 60529)		IP 20
Schutzklasse		Ι
Überspannungskategorie (EN 61010)		II
Verschmutzungsgrad (EN 61010)		2
Nennstrom	A	230 V: Dekontamination: 4,6 Inkubation: 2,4 220 V: Dekontamination: 4,4 Inkubation: 2,3 120 V: Dekontamination: 8,3; Inkubation: 4,6 100 V: Dekontamination: 7,2; Inkubation: 3,9
Leitungsschutzschalter		16 A

Bezeichnung	Einheit	Wert
Nennaufnahme	kW	230 V: Dekontamination: 1,10 Inkubation: 0,56 220 V: Dekontamination: 0,97 Inkubation: 0,51 120 V: Dekontamination: 1,01 Inkubation: 0,55 100 V: Dekontamination: 0,72 Inkubation: 0,39
EMV-Klasse		В

STERI-CYCLE i250

Bezeichnung	Einheit	Wert
Mechanisch		
Außenabmessungen (B x H x T)	mm	780 x 970 x 945
Innenabmessungen (B x H x T)	mm	607 x 670 x 629
Innenraumvolumen, davon Nutzvolumen	1	ca. 255 ca. 162
Einlagebleche (B x T) Anzahl, Lieferumfang Anzahl, maximal Flächenlast, maximal	mm St. St. kg	560 x 500 3 12 10 / Einlageblech
Gesamtlast Gerät, maximal	kg	30
Gewicht, ohne Zubehör	kg	97,5
Thermisch		
Thermische Sicherheitseinrichtungen nach DIN 12880:2007-05		Klasse 3.1 (Temperatur – Wählwächter (TWW) mit Wächterfunktion bei Temperaturüberschreitung)
Umgebungstemperaturbereich	°C	+1834
Umgebungstemperatur gestapelter Geräte	°C	+1828
Temperaturregelbereich	°C	RT + 355
Temperaturabweichung, zeitlich (DIN 12880, Teil 2)	°C	± 0,1
Temperaturabweichung, räumlich (DIN 12880, Teil 2) bei 37 °C $^{*1)}$	°C	± 0,3
Dauer der auto-start-Routine: bis 37 °C Umgebungstemperatur 20 °C	Std.	510
Wärmeabgabe an Umgebung: bei 37 °C während Steri-Run-Dekontamination	kWh/h kWh/h	0,07 0,75
Feuchte		
Wasserqualitäten		Elektrischer Widerstand: 50 kOhmcm bis 1 MOhmcm Leitfähigkeit: 1 bis 20 µS/cm
Füllmengen: Inkubationsbetrieb	1	max. 3 / min. 0,5
Konstantfeuchte bei 37 °C (High-Feuchte Modus) Konstantfeuchte bei 37 °C (Low-Feuchte Modus)	% rF % rF	ca. 93 ca. 90

Bezeichnung	Einheit	Wert
Sonstiges		
Schalldruckpegel (DIN 45 635, Teil 1)	dB(A)	< 50
Relative Feuchte der Umgebung	% rF	max. 80
Höhe des Aufstellungsortes	m NN	max. 2000

^{*1)} Werte ermittelt in Anlehnung an DIN 12880 für Geräte in Standardausführung. Nähere Angaben sind der Kalibrieranweisung zu entnehmen.

STERI-CYCLE i250

Bezeichnung	Einheit	Wert
CO ₂ -Gastechnik		
Gasreinheit	%	min. 99,5 oder med. Qualität
Vordruck	bar	min. 0,8 - max. 1
Mess- und Regelbereich	Vol -%	020
Regelabweichung, zeitlich	Vol -%	± 0,1
CO ₂ -Messzelle		
Genauigkeit bei 37 °C und 5% CO ₂	%CO ₂	± 0,3
O ₂ -Gastechnik		
Gasreinheit	%	min. 99,5 oder med. Qualität
Vordruck	bar	min. 0,8 - max. 1
Mess- und Regelbereich	VoI -%	1 21 bzw. 590
Regelabweichung, zeitlich	VoI -%	± 0,2
O ₂ -Messzelle		
Genauigkeit bei 37 °C und 21% O ₂	%O ₂	± 0,5 (Option: 121% O ₂) ± 2,0 (Option: 590% O ₂)
Elektrisch		
Nennspannung	V	1/N/PE 230 V, AC (± 10%) 1/N/PE 220 V, AC (± 10%) 1/N/PE 120 V, AC (± 10%) 1/N/PE 100 V, AC (± 10%)
Nennfrequenz	Hz	50/60
Schutzart (IEC 60529)		IP 20
Schutzklasse		Ι
Überspannungskategorie (EN 61010)		II
Verschmutzungsgrad (EN 61010)		2
Nennstrom	Α	230 V: Dekontamination: 5,5 Inkubation: 3,3 220 V: Dekontamination: 5,3 Inkubation: 3,2 120 V: Dekontamination: 10,4 Inkubation: 6,3 100 V: Dekontamination: 8,9 Inkubation: 5,3
Leitungsschutzschalter		16 A

Bezeichnung	Einheit	Wert
Nennaufnahme	kW	230 V: Dekontamination: 1,26 Inkubation: 0,76 220 V: Dekontamination: 1,16 Inkubation: 0,69 120 V: Dekontamination: 1,25 Inkubation: 0,75 100 V: Dekontamination: 0,89 Inkubation: 0,53
EMV-Klasse		В

11 Technische Daten STERI-CYCLE i250

Datenkommunikation

Inhalt

- "Struktur der Befehlssequenzen zur Datenkommunikation" auf Seite 12-4
- "Übersicht Allgemeine Parameter (Adressen 0xxx)" auf Seite 12-5
- "Übersicht Inkubator-Parameter (Adressen 2xxx)" auf Seite 12-5
- "Struktur Fehlerspeicher" auf Seite 12-8
- "Struktur Datenlogger" auf Seite 12-11
- "Code-Beispiele Datenlogger" auf Seite 12-15
- "Programm STERI-CYCLE i160 / STERI-CYCLE i250" auf Seite 12-21

USB-Schnittstelle

Die Geräte sind mit einer USB-Schnittstelle ausgerüstet. Die USB-Schnittstelle entspricht dem Standard USB 1.1 / USB 2.0 / USB 3.0 (full speed). Die USB-Schnittstelle wird als virtueller Com Port betrieben. Deshalb kann die Übertragungsgeschwindigkeit der Schnittstelle innerhalb der definierten Baudraten (9.600, 19.200, 38.400, 57.600 Baud) verändert werden. Der Datenaustausch erfolgt über eine festgelegte Struktur von Befehlssequenzen. Die Befehlssequenzen entsprechen dem Aufbauschema der RS 232-Schnittstelle.

Hinweis USB-Anschluss mit virtuellen Com Port einrichten:

Soll zum Datenaustausch zwischen dem PC und dem Inkubator die USB-Schnittstelle verwendet werden, wird mit Hilfe des mitgelieferten Treibers der USB-Anschluss als virtueller Com Port (USB Serial Port) eingerichtet.

Der zugeordnete Com Port wird im Windows-Dialogfenster Gerätemanager/Anschlüsse ermittelt; z. B: USB Serial Port (COM5) und wird dann im Programm **STERI-CYCLE i160 / STERI-CYCLE i250** als Kommunikationsschnittstelle definiert (siehe "Programm STERI-CYCLE i160 / STERI-CYCLE i250" auf Seite 12-21).



Der Treiber ist lauffähig unter den Betriebssystemen: WIN 7, WIN 8, WIN 2000, WIN XP, WIN VISTA.

Treiber USB-Schnittstelle installieren

USB-Kabel an die USB-Schnittstelle (optional) am Schaltkasten des **STERI-CYCLE i160** / **STERI-CYCLE i250** anschließen und mit einem PC verbinden.

Sobald der Hardware-Detektor von Windows den USB-Anschluss erkennt, wird das Dialogfenster ASSISTENT FÜR DAS SUCHEN NEUER HARDWARE geöffnet.

1. Die Option KEINE SOFTWARE SUCHEN wählen.



2. Die Option SOFTWARE VON EINER BESTIMMTEN QUELLE INSTALLIEREN wählen



3. Als Quelle die Daten-CD wählen.



4. Auf der Daten-CD das Unterverzeichnis DRIVER auswählen.

Browse For Folder	?×
Select the folder that contains drivers for your hardware.	
My Documents	^
🗉 😼 PC15	
🖽 🚜 31/2 Floppy (A:)	
🗉 🥯 Festplatte_PC15 (C:)	
🖬 🥝 HERAcell_v01 (D:)	-
🖃 🗁 _Driver	
🗀 amd64	
🗀 i386	
🗀 _Programs	
C DE	
ES ES	
E FR	
🗋 G8	~
To view any subfolders, click a plus sign above.	
ок с	ancel .::

5. Die Installationsroutine installiert den Treiber: EVAL22 Board USB. Nach erfolgreicher Installation wird die Routine mit FERTIGSTELLEN beendet.

Die Übertragungsgeschwindigkeit der Schnittstelle kann innerhalb der definierten Baudraten (9.600, 19.200, 38.400, 57.600 Baud) am Tastbildschirm des **STERI-CYCLE i160 / STERI-CYCLE i250** eingestellt werden (siehe "Einstellungen/Setup" auf Seite 6-17).

Struktur der Befehlssequenzen zur Datenkommunikation

Alle gesendeten und empfangenen Zeichen im Datenaustausch zwischen einem PC und dem Inkubator **Steri-Cycle i160 / Steri-Cycle i250** sind ASCII-Zeichen, die in einem normalen Terminal dargestellt werden können.

Dadurch ist eine einfache Inbetriebnahme, Kontrolle und Programmierung der Kommunikation möglich.

Beschreibung des Protokolls

Zeichencodierung: ASCII-Zeichen, Großbuchstaben sind nicht erlaubt.

Parameter lesen:	
Anfrage:	?:aaaa:bb::cc <cr></cr>
oder:	?:aaaa:bb:XXXX:cc <cr></cr>
Antwort:	!:aaaa:bb:XXXXX:cc <cr></cr>
mit:	aaaa = Adresse des Parameters
	bb =Anzahl der Nutzdaten in diesem Telegramm (00 – ff)
	cc = Checksumme: CRC8-CCITT: $x8 + x2 + x1 + 1 = 0x07$
	ohne cc und <cr></cr>
	XXXX = bb-Bytes Nutzdaten
Beschreibung der Ar	ntwort-Elemente:
aaaa	Adresse des Parameters
ЬЬ	Anzahl der Nutzdaten in diesem Telegramm (00 – ff)
сс	Checksumme: Invertiertes XOR aller Bytes ohne
	Checksumme und <cr></cr>
Beispiel Abfrage der	Softwareversion (50111927):
Abfrage:	?:0001:00::cc <cr></cr>
Antwort:	!:0001:08:50111927:cc <cr></cr>
Parameter schreiben:	
Befehl:	!:aaaa:bb:XXXXX:cc <cr></cr>
Antwort:	!:aaaa:bb::cc <cr></cr>
mit:	aaaa = Adresse des Parameters
	bb =Anzahl der Nutzdaten in diesem Telegramm (00 – ff)
	cc = Checksumme: CRC8-CCITT: $x8 + x2 + x1 + 1 = 0x07$
	ohne cc und <cr></cr>
	XXXX = bb-Bytes Nutzdaten
Antworten mit einer	Fehlermeldung:
Antwort:	!:aaaa:bb:XX:cc <cr></cr>
Beschreibung	g der Antwort-Elemente:
--------------	--
aaaa	Adresse des Parameters,
ЬЬ	Anzahl der Nutzdaten (immer 02)
сс	Checksumme: CRC8-CCITT: x8 + x2 + x1 + 1 = 0x07
	ohne cc und <cr></cr>
	XX = 2 Bytes Fehlermeldung (siehe Tabelle unten)

Beispiel unbekannter Befehl:				
Anfrage:	?:0005:00::cc <cr></cr>			
Antwort	!:0005:02:?1:cc <cr></cr>			

Bedeutung der beiden Bytes in der Fehlermeldung:

Fehlermeldung	Beschreibung
?0	Fehler in der Telegrammstruktur oder Checksumme
?1	Unbekannter Befehl oder unbekannter Parameter
?2	Interner Speicherfehler
?3	Datenfehler (Wert nicht innerhalb seiner Grenzen)

Übersicht Allgemeine Parameter (Adressen 0xxx)

Allgemeine Parameter sind Systemwerte, wie Datum, Uhrzeit und die Versionsnummer der Hauptplatine.

Allgemeine Parameter lesen

Adresse	Beschreibung	Bemerkung
0001	Versionsnummer Mainboard	8 Stellen
0010	Ausgabe Datum und Uhrzeit [Stunden:Minuten:Sekunden]; [Tag:Monat:Jahr]	17 Bytes / Dezimalwert im Format xx:xx:xx;xx:xx
0011	Datum [Tag:Monat:Jahr]	8 Bytes / Dezimalwert im Format xx:xx:xx
0012	Uhrzeit [Stunden:Minuten:Sekunden]	8 Bytes / Dezimalwert im Format xx:xx:xx

Übersicht Inkubator-Parameter (Adressen 2xxx)

Inkubator-Parameter sind unterteilt in:

- die Parameter (Basic) der drei Regelkreise Temperatur, CO₂ und O₂,
- die Parameter (interne Funktionen) der Betriebsfunktionen und Datenaufzeichnung.

Parameter (Basic) lesen

Adresse	Beschreibung	Bemerkung
2000	Gerätestatus ^{*1)} (Fehler-) Status der Regelkreise Temperatur, CO ₂ , O ₂ , rH, Ref-Temp.	33 Bytes / Hexadezimalwert im Format xxxxxxxx;xxxx; ;xxxx;xxxx;xxxx
2010	Soll-, Ist- und Referenztemperatur ^{*2})	23 Bytes / Dezimalwert im Format +xxx.xx;+xxx.xx;+xxx.xx
2020	Soll- und Ist-CO ₂ -Gehalt ^{*2)}	15 Bytes / Dezimalwert im Format +xxx.xx;+xxx.xx
2030	Soll- und Ist-O ₂ -Gehalt ^{*2)}	15 Bytes / Dezimalwert im Format +xxx.xx;+xxx.xx
204a	Ist-Wasserstand (100% oder 0%)	7 Bytes / Dezimalwert im Format +xxx.xx
204b	Anzeige Low-Feuchte (1 aktiv, 0 nicht aktiv)	2 Bytes / Hexadezimal im Format xx

^{*1)} Beispiel Gerätestatus und (Fehler-) Status Regelkreise (Details siehe Fehlermeldungen)

^{*2)} Alle Werte besitzen 2 Nachkommastellen

Parameter	(interne	Funktionen)) lesen
-----------	----------	-------------	---------

Adresse	Beschreibung	Bemerkung
2100	Status Ablauf ^{*1)} und Restlaufzeit [Stunden:Minuten] Desinfektion, sowie Datum und Uhrzeit des letzten Starts	25 Bytes / Dezimalwert im Format xx;+xxx:xx;xx.xx;xx:xx
2105	Status Ablauf ^{*1)} , aktueller CO ₂ Offset + Wartezeit [Minuten:Sekunden] auto-start, sowie Datum und Uhrzeit des letzten Starts	25 Bytes / Dezimalwert im Format xx;xx.x;+xxx:xx;xx.xx;xx:xx
2140	Status Gasflaschenumschalter CO ₂ lesen ^{*3)}	2 Bytes / Hexadezimal im Format xx
2141	Status Gasflaschenumschalter O ₂ lesen ^{*3)}	2 Bytes / Hexadezimal im Format xx
2300	Fehlerspeicher auslesen (aktuelle Fehler) ^{*4)}	Bis zu 241 Bytes / Hexadezimalwert Format siehe eigenes Kapitel
2301	Fehlerspeicher auslesen (ältere Fehler) ^{*4)}	Bis zu 241 Bytes / Hexadezimalwert Format siehe eigenes Kapitel
2400	Abfrage (Start) der im Datenlogger gespeicherten Daten ^{*5)}	Bis zu 224 Bytes / Hexadezimalwert Format siehe eigenes Kapitel
2401	Abfrage der weiteren im Datenlogger gespeicherten Daten ^{*6)}	224 Bytes / Hexadezimalwert Format siehe eigenes Kapitel
2402	(Wiederholungs-) Abfrage der letzten Datenloggerabfrage ^{*7)}	224 Bytes / Hexadezimalwert Format siehe eigenes Kapitel
2410	Schreibzyklus Datenlogger in Stunden:Minuten:Sekunden auslesen	8 Bytes / Dezimalwert im Format xx:xx:xx

*1) Siehe Tabelle Anmerkungen Status Ablauf Desinfektion und auto-start.

*2) Jeweils 2 Bytes pro Ebene.

*3) Flasche A aktiv (0x01), Flasche B aktiv (0x02), Druck Flasche A OK (0x10), Druck Flasche B OK (0x20).

*4) Weitere Informationen über den Fehlerspeicher in Kapitel 13.5.

*5) Lesezeiger auf den ersten Eintrag setzen, max. 7 Einträge lesen.

*6) Die nächsten 7 Einträge senden. Lesezeiger automatisch auf den nächst neueren Eintrag setzen, max. 7 Einträge lesen.

*7) Einträge des letzten Telegramms nochmals senden. Kann genutzt werden, wenn die Kommunikation gestört war.

Bit	Desinfektion	auto-start
0x00	Steri-run nicht aktiviert	auto-start nicht aktiviert
0x01	Initialisierung	Initialisierung
0x02	Türöffnungszeit abwarten	Türöffnungszeit abwarten
0x03	Türschließen abwarten	Türschließen abwarten
0x04	Start	Start
0x05	Heizen	Heizen
0x06	Halten	Gegenspannungsableich durchführen
0x07	Kondensation	Wartezeit 1
0x08	Kühlen	Toleranzband setzen
0x09	Trocknen	Stabile Feuchte aufbauen
0x0A	Warten auf Freigabe	Gegenspannungsableich durchführen
0x0B	Abbruch	Wartezeit 2
0x0C	-	Offset ermitteln
0x0D	-	Offset auslesen, prüfen
0x0E	-	Freigabe
0x0F	-	Abbruch

Anmerkung zu *3) Status Ablauf Desinfektion und auto-start:

Struktur Fehlerspeicher

Der Fehlerspeicher enthält 22 Fehlermeldungen. Die Abfrage wird mit jeweils 22 Datensätzen, die mit einem Doppelpunkt als Trennzeichen getrennt werden, beantwortet und kann mit den folgenden Befehlen abgefragt werden:

Abfrage:	?:2300:00::cc <cr> Lesen der letzten 11 Einträge des Fehlerspeichers.</cr>
Abfrage:	?:2301:00::cc <cr> Lesen der ersten 11 Einträge des Fehlerspeichers.</cr>

Diese Datensätze bestehen aus 11 Byte und werden vor der Übertragung in 21 ASCII-Zeichen verschlüsselt. Z.B. werden aus dem Byte 0x23 die ASCII-Zeichen 0x32 ("2") und 0x33 ("3").

- Byte 1 besteht aus 1 Zeichen,
- Bytes 2 11 bestehen aus 2 Zeichen.

Somit besteht eine Antwort aus 1+(10 x 2) = 21 Daten-Bytes plus Trennzeichen. In einem Datensatz wird immer das Datum, die Uhrzeit, der fehlerhafte Regelkreis, der Gerätestatus und die Fehlermeldung übergeben.

Antwortbeispiel:	
!:2300:fb:10b01060	f372280000002:20b01060f38100001:80
Erster Datensatz:	!:2300:fb:10b01060f372280000002: (mit 21 Bytes)
Zweiter Datensatz:	20b01060f38100001:80 (Beginn des zweiten Datensatzes nach 01060 Bytes des ersten Datensatzes und des Trennzeichens [1 Byte])

Schema Datensatzstruktur Fehlerspeicher:

	00	0b	01	06	0f	37	22	8000	0002
Typ fehlerhafter Regelkreis									
0x00 = Fehler im									
Regelkreis Temperatur	00								
Datum / Tag									
0x0b = der 11. des Monats		0b							
Datum / Monat									
0x01 = Januar			01						
Datum / Jahr									
$0 \times 06 = 2006$				06					
Uhrzeit / Stunden					-				
0x0f = 15 Uhr					0f				
Uhrzeit / Minuten									
0x37 = 55 Minuten						37			
Uhrzeit / Sekunden									
0x22 = 34 Sekunden							22		
Gerätestatus									
0x8000 = Gerätefehler vor-									
handen								8000	
Fehlermeldung									
0x0002 = lstwert über									0002

In diesem Datensatz werden die folgenden Informationen übertragen:

- Erstellt am 11. Januar 2006 um 15:55:34 Uhr.
- Ein Gerätefehler liegt an und die Ist-Temperatur ist zu hoch.

Übersicht der möglichen Fehlermeldungen in Hex-Codierung

Hex-Code	Beschreibung / Typ
0x00	Regelkreis Temperatur
0x01	Regelkreis CO ₂
0x02	Regelkreis O ₂
0x07	Wasserstand
0x08	Allgemeiner Gerätestatus

Übersicht der möglichen Fehlermeldungen in Bit-Codierung Allgemeiner Gerätestatus, Regelkreis Temperatur und CO₂:

Bit	Allgemeiner Gerätestatus
0x0002	Gerätetür zu lange offen
0x0004	Display kommuniziert nicht
0x0008	Parameter Mainboard unplausibel (EEPROM defekt)
0x0010	Datenlogger defekt (Gerät ist weiter funktionsfähig)
0x0020	Fehler in der Desinfektion / Steri-run
0x0040	Keine Stromversorgung während Steri-run
0x0080	Fehler im auto-start
0x0100	ADC Test fehlgeschlagen
0x0400	Lüfterfehler
0x1000	IR Sensor getauscht (Info)
0x2000	auto-start aktiv (Info)
0x4000	Desinfektion aktiv (Info)
0x8000	Gerätefehler vorhanden (Info)
Bit	Fehlerstatus Regelkreis Temperatur
0x0001	Sensor-/ Fühlerbruch
0x0002	Istwert über
0x0004	Istwert unter
0x0008	Istwert nicht plausibel
0x0010	Kalibrierwerte zu groß/klein

Bit	Fehlerstatus Regelkreis CO ₂
0x0001	Sensor-/ Fühlerbruch
0x0002	Istwert über

Bit	Fehlerstatus Regelkreis CO ₂
0x0004	Istwert unter
0x0010	Kalibrierwerte zu groß/klein
0x0020	Störung Kommunikation (zum Sensor)
0x0040	Störung Kommunikation (zum Gasflaschenumschalter)
0x0080	Kein Gas vorhanden, Flasche A und B leer
0x0200	Gasflasche A leer
0x0400	Gasflasche B leer

Regelkreis O₂ und Wasserstand:

Bit	Fehlerstatus Regelkreis O ₂
0x0001	Sensor-/ Fühlerbruch
0x0002	Istwert über
0x0004	Istwert unter
0x0020	Störung Kommunikation (zum Sensor)
0x0040	Gasflaschenumschalter kommuniziert nicht
0x0080	Kein Gas vorhanden, Flasche A und B leer
0x0200	Gasflasche A leer
0x0400	Gasflasche B leer
Bit	Fehlerstatus Wasserstand
0x0001	Kein Wasser

Struktur Datenlogger

Der Datenlogger speichert bis zu 10.000 Einträge. Abhängig von der Einstellung des Speicherzyklus (in Sekundenabschnitten) kann - z. B. bei einem Wert von 10.000 s (Defaultwert) - der Speicherzeitraum die Ereignisse von ca. 5 Tagen aufzeichnen.

Im Datenlogger werden folgende Informationen gespeichert:

- Wichtige Benutzeraktionen, Systemereignisse und Fehlermeldungen.
- Messdaten der drei Regelkreise während des Inkubationsbetriebes.

Der Datenlogger kann mit den folgenden Befehlen abgefragt werden:

?:2400:00::cc <cr></cr>
Setzen des Lesezeigers des Datenloggers auf den ältesten Eintrag und Ausgabe der
ersten Datensätze.
?:2401:00::cc <cr></cr>
Ausgabe der folgenden Datensätze, der Lesezeiger wandert automatisch
schrittweise von den älteren zu den aktuellen Einträgen.

Abfrage:	?:2402:00::cc <cr></cr>		
	Erneute Ausgabe der zuletzt gelesenen Daten, der Lesezeiger wird bei diesem		
	Befehl nicht verschoben. Mit diesem Befehl kann ein Datenverlust nach		
	einem Kommunikationsfehler verhindert werden.		
Die Abfragebefehl	e werden jeweils mit bis zu 7 Datensätzen, die ohne ein Trennzeichen aufeinander		
folgen, beantworte	et. Diese Datensätze bestehen aus 16 Bytes und werden vor der Übertragung in 32		
ASCII-Zeichen ve	rschlüsselt.		
So werden z.B. aus	dem Byte 0x23 die ASCII-Zeichen:		
0x32 ("2") und 0x33 ("3").			
Folglich besteht eine Antwort aus bis zu 7 x 16 = 112 Bytes und somit 224 ASCCII-Zeichen.			
In einem Datensatz wird immer das Datum und die Uhrzeit (ohne Sekunden), der Gerätestatus und			
der Typ des Datenloggereintrags übergeben (Byte 0-7, bzw. ASCII-Zeichen 0-15).			
Je nach Eintrag können weiterhin die aktuellen Ist- oder Sollwerte der Regelkreise oder sonstige			
Parameter eingetragen werden (Byte 8-15, bzw. ASCII-Zeichen 16-31).			
Antwortbeispiel:			
1.2400:e0.010b01060f3700000177002800d40000110b01060f3800000172003200d20352			

!:2400:e0:010b01060f3700000177002800d4000011
(bestehend aus 32 Byte ASCII-Zeichen)
0b01060f3800000172003200d20352:80
(Beginn des zweiten Datensatzes nach 32 Bytes des ersten Datensatzes)

Schema Datensatzstruktur Datenlogger:

Typ Datenloggereintrag	01	0b	01	06	Of	37	0000	0177	0028	00d4	0000
0x01 = Standardeintrag der alle 60 s gespeichert wird	01										
Datum / Tag 0x0b = der 11. des Monats		0b									
Datum / Monat 0x01 = Januar Datum / Jahr			01								
0x06 = 2006 Uhrzeit / Stunden 0x0f = 15 Uhr				06	Of						
0x37 = 55 Minuten Gerätestatus						37					
Typabhängig, hier aktuelle Temperatur in 1/10 °Celsius 0x0177 = 37,5°C							0000	0177			
Typabhängig, hier aktuelle CO2-Gehalt in 1/10% 0x0028 = 4,0% CO ₂									0028		
Typabhängig, hier aktuelle O₂-Gehalt in 1/10%											
$0x00d4 = 21,2\% O_2$										00d4	
Reserviert für rH in 1/10% immer 0x0000 = 0%											0000

In diesem Datensatz werden die folgenden Informationen übertragen:

- Erstellt am 11. Januar 2006 um 15:55 Uhr.
- Der Gerätestatus meldet keine Besonderheiten.
- Temperatur beträgt 37,5 °C.
- Gaskonzentration 4,0 % CO₂, 21,2 % O₂.

Hinweis Code-Beispiel:

Ein Code-Beispiel finden Sie im Anschluss an dieses Kapitel.

Übersicht der möglichen Ereigniseinträge in Bit-Codierung Übersicht Ereigniseinträge Teil I:

Code	Ereignis	Informationen (Byte 8-15)				
0x01	Sollwerte von allen Regelkreisen (periodisch in Minutenzyklen)	Aktuelle Werte von Temperatur, CO ₂ , O ₂ und rH				
0x02	Sollwertänderung (am Beginn eines neuen Abschnittes)	Sollwert von Temperatur, CO ₂ , O ₂ und rH				
0x10	Sollwert Temperatur ändern	Sollwert von Temperatur, CO ₂ , O ₂ und rH				
0x11	Sollwert CO _{2 ändern}	Sollwert von Temperatur, CO ₂ , O ₂ und rH				
0x12	Sollwert O _{2 ändern}	Sollwert von Temperatur, CO ₂ , O ₂ und rH				
0x20	Neuer Fehler Temperatur	Status / Fehlerregistrierung bei Temperatur, CO ₂ , O ₂ und rH				
0x21	Neuer Fehler CO ₂	Status / Fehlerregistrierung bei Temperatur, CO ₂ , O ₂ und rH				
0x22	Neuer Fehler O ₂	Status / Fehlerregistrierung bei Temperatur, CO ₂ , O ₂ und rH				
0x2F	Neuer Fehler System	Status / Fehlerregistrierung bei Temperatur, CO ₂ , O ₂ und rH				
0x30	Netzreset	Sollwert von Temperatur, CO ₂ , O ₂ und rH				
0x31	Deckel offen	Aktuelle Istwerte von Temperatur, CO ₂ , O ₂ und rH				
0x32	Tür geschlossen	Aktuelle Istwerte von Temperatur, CO ₂ , O ₂ und rH				
0x40	Kundenkalibration Temperatur	Kalibrationsebene (2 Bytes), alte Temperatur, neue Temperatur (2 Bytes jede)				
0x41	Kundenkalibration CO ₂	Kalibrationsebene (2 Bytes), alter CO ₂ -Wert, neuer CO ₂ -Wert (2 Bytes jede)				
0x42	Kundenkalibration O ₂	Kalibrationsebene (2 Bytes), alter O ₂ -Wert, neuer O ₂ -Wert (2 Bytes jede)				
0x50	Start auto-start	Status / Fehlerregistrierung bei Temperatur, CO ₂ , O ₂ und rH				
0x51	auto-start erfolgreich beendet	Aktuelle Istwerte von Temperatur, CO ₂ , O ₂ und rH				

Code	Ereignis	Informationen (Byte 8-15)
0x52	auto-start beendet mit Fehler	Status / Fehlerregistrierung bei Temperatur, CO ₂ , O ₂ und rH
0x53	auto-start manuell gestoppt	Status / Fehlerregistrierung bei Temperatur, CO ₂ , O ₂ und rH
0x60	Start Steri-run	Status / Fehlerregistrierung bei Temperatur, CO ₂ , O ₂ und rH
0x61	Steri-run erfolgreich beendet	Aktuelle Istwerte von Temperatur, CO ₂ , O ₂ und rH
0x62	Steri-run beendet mit Fehler	Status / Fehlerregistrierung bei Temperatur, CO ₂ , O ₂ und rH
0x63	Steri-run manuell gestoppt	Status / Fehlerregistrierung bei Temperatur, CO ₂ , O ₂ und rH
0x70	Gas Monitor Flasche A leer	Status Gasmonitoring (2 Bytes), 4 Bytes leer
0x71	Gas Monitor Flasche B leer	Status Gasmonitoring (2 Bytes), 4 Bytes leer
0x72	Gas Monitor manuelle Umschaltung	Status Gasmonitoring (2 Bytes), 4 Bytes leer
0x90	Start Low Feuchte	Aktuelle Istwerte von Temperatur, CO ₂ , O ₂ und rH
0x91	Stop Low Feuchte	Aktuelle Istwerte von Temperatur, CO ₂ , O ₂ und rH
0xe0	Löschen des Datenloggers	Aktuelle Istwerte von Temperatur, CO ₂ , O ₂ und rH
0xff	Letzter Eintrag des Datenloggers	Keine Information, auch keine Information zu Datum, Zeit und Status

Übersicht Ereigniseinträge Teil II:

Code-Beispiele Datenlogger

Ein Eintrag im Datenlogger hat eine Größe von 16 Bytes und ist wie folgt aufgebaut:

- 1. Byte: gibt das Ereignis an (z.B. Tür geöffnet 0x31, Messwerteintrag 0x01)
- 2. Byte: Tag des Eintrags
- 3. Byte: Monat
- 4. Byte: Jahr
- 5. Byte: Stunde
- 6. Byte: Minuten
- 7. und 8. Byte: Gerätestatus
- 9. bis 16. Byte: verschiedenste Daten zum Ereignis

Funktionen zur Abfrage Datenlogger

Im folgenden Code-Beispiel zum Auslesen des Datenloggers werden sechs Funktionen benutzt:

- ahex
 - // wandelt das empfangene ASCII-Zeichen in ein Hexadezimalzahl um,
- send_telegramm
 // sendet eine Anfrage an den Datenlogger,
 oot telegramm
- get_telegramm
 // empfängt eine Antwort vom Datenlogger,
- time_2_str // erstellt aus einem Hex-Wert ASCII-Zeichen im Uhrzeitformat,
- read_datalogger
 // bearbeitet die empfangenen Daten und schreibt diese in eine Datei.

Code-Beispiel zur Abfrage Datenlogger

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;
// Telegramm zusammenkopieren
    strncpy (&string[0], ":xxxx:00::00\r", 14);
// 4-stellige Adresse einfügen
    strncpy (&string[2], p, 4);
// Checksumme berechnen: invertiertes XOR aller Bytes
// ohne Checksumme und <CR>
    for (i = 0; i < 11; i + +)
        bcc = (bcc^string[i]);
// Checksumme kopieren
    string[11] = hexa(bcc/16);
    string[12] = hexa(bcc%16);
```

```
// Telegramm senden
ComWrt (COM_NR, string, 14);
return;
}
```

get_telegramm

```
int get_telegramm(char *p)
{
    int reading_count = 0;
// zeichenweise Auslesen des Telegramms
    do
        ComRd(COM_NR, &p[reading_count], 1);
// bis zum Erhalt von <CR>
    while ((p[reading_count++] != '\r'));
// Rückgabe = Anzahl der empfangenen
    Zeichenreturn (reading_count);
```

time_2_str

```
char time_2_str (int z, char * b)
{
     char i;
// zwei Ziffern ausgeben
     for (i = 1; i >= 0; i--){
//Wert berechnen
        b[i] = z%10+0x30;
// Vorgabewert verkleinern
        z = z/10;
    }
    return (2);
    }
```

num_2_string

```
char num_2_str (int z, char * b)
{
// Zahl mit einer Nachkommastelle
    char a[12];
    char i, l;
    int rest = 0;
    l = 0;
// Negative Zahl?
    if (z < 0) {
// Vorzeichen setzen
        b[0] = '-';l = 1;
// Wert umrechnen
        z = 0xffffffff-z+1;
    }
// Nachkommastelle speichern</pre>
```

```
rest = z \% 10;
// Nachkommastelle abschneiden
    z = z / 10;
// Zahl vor dem Komma berechnen und kopieren
    for (i = 0; i < 12; i++){
// Wert berechnen
    a[i] = z\%10 + 0x30;
// Vorgabewert verkleinern
    z = z/10;
// Zahl komplett kopiert?
    if (z == 0) break;
    ł
    for (; i \ge 0; i--)b[l++] = a[i];
// Zahl nach dem Komma berechnen und kopieren
    b[l++] = ',';
// Wert berechnen
    b[l++] = rest\%10+0x30;
    return (l);
    ł
```

read_datalogger

```
int read_datalogger ()
#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,
// schreiben der Titelzeile in der Datei
WriteFile (FileHandle, "Date; Time; Comment; Temp Act.; CO2 Act.; O2
Act.;rH Act.;Temp Set;CO2 Set;O2 Set;rH Set;\n", 85);
// Endlosschleife
while (1)
// Datenlogger auf Anfang setzen und lesen
    if (!GetTele) {
    send_telegramm ("2400");
    }
    else{
// weitere Datensätze lesen
    send_telegramm ("2401");
    len = get_telegramm (buffer);
// kein Telegramm erhalten
```

```
if (!len) \{
        GetError ++;
// erneute Anfrage
    send_telegramm ("2402");
    len = get_telegramm (buffer);
// wieder kein Telegramm erhalten
    if (!len) return 1;
    }
// Telegrammzaehler erhöhen
    GetTele ++;
// Länge der gesendeten Nutzdaten
    len = (ahex(buffer[7]) * 0x10 + ahex(buffer[8])) / 2;
// Wandlung ASCII String in einen nutzbaren Zahlenstring
    for (i = 0; i < (string); i++)
    zahlenstring [i] = (ahex(buffer[10 + (2*i)]) * 0x10 +
    ahex(buffer[11 + (2*i)]));
// Berechnung der gesendeten Datenpakete
    data = ((len) / SIZE_DATA2);
// Auswertung aller Datenpakete
    for (i = 0; i < data; i++)9
    len = 0;
// Uhrzeit und Datum in Datei schreiben
    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:
//zyklische Einträge prüfen auf Gerätefehler
    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{
// alle Gerätefehler abfragen (siehe "Übersicht der möglichen Ereigniseinträge in Bit-Codierung
```

```
Übersicht Ereigniseinträge Teil I:" auf Seite 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;
                     }
                }
// jetzt restliche Gerätefehler abfragen
11
//
11
11
//
//und schließlich zyklischen Istwerteintrag ohne Gerätefehler
//abfragen
else{
            str_cpy (&string[string], ",ok;", 3);
            string += 3;
        }
    1
// Istwerte aus Zahlenstring in Excelstring kopieren
    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 ++] = ';';
// ab hier Sollwerte eintragen
    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;
```

```
// ab hier Abfrage der restlichen Events
    case EVENT_FORMAT_DATALOG:
        WriteFile (FileHandle, excelstring, len);
        WriteFile (FileHandle, "Data logger erased;\n",20);
        break;
    case EVENT_POWER_ON:
// Aktualisierung der Sollwerte
        SetTemp = zahlenstring [8+i*SIZE_DATA2]*0x100+
        zahlenstring[9+i*SIZE_DATA2];
        SollCO2 = zahlenstring[10+i*SIZE_DATA2]*0x100+
        zahlenstring[11+i*SIZE_DATA2];
        SollO2 = 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 ...
//hier alle Events abfragen (siehe "Übersicht der möglichen Ereigniseinträge in Bit-Codierung
Übersicht Ereigniseinträge Teil I:" auf Seite 12-14)
// Abbruch 0xFF kennzeichnet das Ende des Datenloggers
    case 0xFF:
    WriteFile (FileHandle, "End;\n",5);
    }
}
return 0;
```

Programm STERI-CYCLE i160 / STERI-CYCLE i250

Das Programm stellt eine Benutzeroberfläche (nur mit englischen Menübezeichnungen) zur Abwicklung der Datenkommunikation zwischen dem Gerät und einem angeschlossenen PC zur Verfügung.

1	Thermo Fisher CO2 Incubators
	Thermo Fisher CO2 Incubators
	It is strongly recommended that you exit all programs before running this installer. Applications that run in the background, such as virus-scanning utilities, might cause the installer to take longer than average to complete.
	<< Back Next >> Cancel

Das Programm dient zum:

- Auslesen und Archivieren von Fehlermeldungen (Error Logger). Die Datensätze werden in dem Metaformat *.CSV gespeichert.
- Auslesen und Archivieren von Ereigniseinträgen (data Logger). Die Datensätze werden in dem Metaformat *.CSV gespeichert.
- Erstellen einer Service-Datei (Servicefile) zum Versand an den Technischen Service von Thermo Fisher Scientific. Anhand der Informationen der Service-Datei kann eine systematische Fehlerermittlung durchgeführt werden. Die Datensätze werden in dem proprietären Format *.SRF gespeichert:

STERI-CYCLE i160 & STERI-CYCLE i250 installieren

- 1. Installationsroutine starten:
- Auf der Daten-CD im Unterverzeichnis PROGRAMS die Datei SETUP.EXE mit Doppelklick aufrufen.

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

2. Installationsverzeichnis für das Programm bestimmen.

Select the primary installation directory.		
All software will be installed in the followin different location, click the Browse buttor	g locations. To install softw and select another directo	vare into a rry.
Target directory for application C:\Program Files (x86)\TFS\		Browse
Target directory for National Instrument	s software	
C:\Program Files (x86)\National Instrum	ients\	Browse

- 3. In der Abfolge der vorgegebenen Installationsschritte:
 - die Lizenzvereinbarung bestätigen,
 - den Installationsumfang bestätigen,
 - nach Ausgabe der Vollständigkeitsmeldung zur Installation die Installationsoberfläche schließen und den Computer neu starten.

STERI-CYCLE i160 & STERI-CYCLE i250 bedienen

Struktur der Benutzermenüs:

Die Benutzeroberfläche ist gegliedert in zwei Hauptmenüs:

- MAIN mit beiden Funktionselementen:
 - Ausgabe der Programmversion: FIRMWARE VERSION
 - Schalter zum Beenden des Programmes: QUIT

Thermo Fisher CO2 Incubators COM: NoCOM	x
MAIN	GENERAL
50135297 V2.7 20.05 2014 Device Name no com.	ThermoFisher SCIENTIFIC
Date: no com Firmware version: no com	Que

- GENERAL mit den Untermenüs:
 - PRESETTING zur Einstellung der Übertragungsgeschwindigkeit und zur Auswahl des seriellen Anschlusses,
 - TEST COM zum Testen der Kommunikationsverbindung PC zum Inkubator,
 - DATE & TIME zur Einrichtung des Datums und der Zeit auf die gewünschte Zeitzone,
 - ERROR LOGGER zum Auslesen der Fehlermeldungen,
 - DATA LOGGER zum Auslesen der Ereigniseinträge,
 - SERVICEFILE zum Auslesen der Fehlerinformationen und zur Erstellungen einer Service-Datei,
 - PASSWORD sperrt den Zugriff auf die Geräteparameter des Inkubators.

Funktion der Benutzermenüs:

PRESETTING

Das Untermenü PRESETTING ermöglicht die Einstellung der Übertragungsgeschwindigkeit und die Auswahl des seriellen Anschlusses.

Thermo Fisher CO2 Incubators COM: NoCOM						
	MAIN			GENERAL		
PRESETTING	TEST COM	DATE & TIME	LOGGER	SERVICEFILE	PASSWORD	
PRESETTING Serial Baudrate 57 600 V chang save setting	MAIN TEST COM	DATE & TIME	LOGGER	SERVICEFILE	PASSWORD	

- 1. Übertragungsgeschwindigkeit auswählen im Bereich von 9600 115200 Baud.
- 2. Seriellen Anschluss des PC auswählen. Ist der USB-Treiber installiert, kann der dem USB-Anschluss zugeordnete (virtuelle) Com Port ausgewählt werden (siehe "USB-Schnittstelle" auf Seite 12-1).
- 3. Einstellungen übernehmen:
- Taste CHANGE drücken.
- 4. Einstellungen abspeichern (in einer Ini-Datei):
- Taste SAVE TO FILE drücken.

Hinweis Übertragungsgeschwindigkeiten:

Die Einstellungen der Übertragungsgeschwindigkeiten im Benutzermenü PRESETTING und am Gerät müssen identisch sein!

TEST COM

Das Untermenü TEST COM dient zum Testen der Kommunikationsverbindung mit den im Untermenü PRESETTING definierten Einstellungen.

MAIN			GENERAL	
PRESETTING TEST COM	DATE & TIME	LOGGER	SERVICEFILE	PASSWORD
command (!) / request (?) !? address (e.g. 2010) 2010 attachment (e.g. +037.00;) string sent				send
string received				1

- 1. Beispielabfrage der aktuell messbaren Temperaturwerte des Inkubators:
- Abfrage: ? (Voreinstellung, nicht veränderbar)
- Adresse: 2010 (Adresse Temperaturwerte: Sollwert, Istwert, Referenzwert)
- 2. Abfrage an den Inkubator absenden:
- Taste SEND drücken.
- Sendet der Inkubator einen Antwortstring zurück, steht die Kommunikationsverbindung zum Inkubator.
- Ist keine Verbindung möglich, wird ein Fehlerdialog angezeigt:



- 3. Fehlerdialog schließen:
- Taste OK drücken.

DATE & TIME

Das Untermenü DATE & TIME ermöglicht die Anpassung des Datums und der Zeit auf die gewünschte Zeitzone.

MAIN			GENERAL	
PRESETTING TEST COM	DATE & TIME	LOGGER	SERVICEFILE	PASSWORD
device date: 03.07.14 set (6d.mm.yy) device time: 10.30.44 set (hh.mm.sa) system time: 07.07.14 10.55.09 set (dd.mm.yy hh.mm.sa)				

- 1. Die Eingaben in den beiden Textfeldern müssen im Format TT.MM.JJ (Tag, Monat, Jahr) erfolgen.
- 2. Eingabe übernehmen:
- Taste SET drücken.

ERROR LOGGER

Das Untermenü ERROR LOGGER dient zum Einlesen der Fehlermeldungen in das Textfeld der Benutzeroberfläche.

Thermo Fisher CO2 Incubators O	COM: NoCOM				
	MAIN			GENERAL	
PRESETTING	TEST COM	DATE & TIM	E LOGGER	SERVICEFILE	PASSWORD
DATA LOGGER	ERROR LOGGER	1			
	date	time	Control loop	L	oop message
read Amount data 0 Amount telegrams 0 Amount erons 0					: save to file

Die Datensätze können im Metaformat *.CSV abgespeichert werden.

- Datensätze als Datei abspeichern:
- Taste SAVE TO FILE drücken.

DATA LOGGER

Das Untermenü DATA LOGGER dient zum Einlesen der Ereigniseinträge in das Textfeld der Benutzeroberfläche.

Die Datensätze werden im Metaformat *.CSV abgespeichert.

- 1. Datensätze einlesen:
- Taste READ drücken.
- 2. Der Verlauf der Datenübertragung wird in den drei Textfeldern angezeigt:
- AMOUNT DATA: Anzahl der übertragenen Datensätze insgesamt
- AMOUNT TELEGRAMS: Anzahl der übertragenen Telegramme.
- AMOUNT ERRORS: Anzahl der übertragenen Fehlermeldungen.

Hinweis Dauer der Datenübertragung:

Da der Datenlogger bis zu 10.000 Datensätze beinhalten kann, kann die Datenübertragung auf einen PC einige Zeit in Anspruch nehmen.

SERVICEFILE

Das Untermenü SERVICEFILE dient zum Einlesen der Fehlerinformationen des Inkubators und zur Erstellung einer Service-Datei, abgespeichert im proprietären Format *.srf. Die Service-Datei wird zur Fehleranalyse an den Technischen Service von Thermo Fisher Scientific weitergeleitet.

	MAIN				GENERAL	
PRESETTING	TEST COM	DATE & TIME	LOGGI	ER S	SERVICEFILE	PASSWOR
	Öffnen				×	
	Directory					
	History: 10.1					
	Suchen in:	SDisk (C:)	•	+ 🗈 💣 📰 🔻		
	æ	Name		Änderungsdatum	Typ 🔺	
		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	Doku		11.06.2014 13:54	Dateio E	
mount data	Desktop	download		24.01.2014 11:54	Dateio	
)		DPSEDriver		14.11.2013 14:25	Dateio	
	Bibliotheken	drivers		08.01.2014 07:13	Dateio	
mount telegrams		FILES		22.03.2012 13:17	Dateio	
)	1	📕 Hotfix		10.02.2012 00:14	Dateio	
	Computer	퉬 Intel		17.01.2011 08:44	Dateio	
		📕 Kpcms		16.03.2012 08:01	Dateio	
	Netzwerk	🔁 PDM-DELSB		28.09.2012 07:20	Dateio	
	HOLEWOIK	Perfl ons		14 07 2009 05-20	Datein	
		Dateiname: service.srf		_	ОК	
		Dateityp: (*.srf)		•	Abbrechen	

- 1. Service-Datei erstellen:
- Taste CREATE SERVICEFILE drücken.
- Im Windows Dialogfenster den Dateinamen und das Verzeichnis zum Speichern festlegen.
- 2. Speichervorgang starten:
- Taste OK drücken.

Hinweis Dauer der Erstellung:

Die Zusammenstellung der Geräteinformationen und das Erstellen der Service-Datei kann einige Zeit in Anspruch nehmen.

PASSWORD

Das Untermenü PASSWORD steht nur dem Servicepersonal von Thermo Fisher Scientific zur Verfügung.

Thermo Fisher CO2 Incubators COM: NoCOM				
MAIN			GENERAL	
PRESETTING TEST COM	DATE & TIME	LOGGER	SERVICEFILE	PASSWORD
	PASSWORD	ENTER		

13-

Gerätebuch

Gerätetyp:			Bestell-Nr.:		
Fabrik-Nummer:			Service-Nummer:		
Aufstellungsort:			Vermerke des Betreibers:		
Durchgeführte Arbeit	en:	Bemerkunger	n:	Datum	Unterschrift

Bitte die durchgeführten Arbeiten hier eintragen:

13 Gerätebuch

Kontaktdaten Thermo Scientific

Übersicht der internationalen Thermo Fisher Vertriebsorganisationen Postanschrift Deutschland: Thermo Electron LED GmbH Robert-Bosch-Straße 1 D - 63505 Langenselbold

Anfragen aus Deutschland:

Telefon Vertrieb 0800 1 536376 Telefon Service 0800 1 112110 Fax Vertrieb/Service 0800 1 112114 E-Mail info.labequipment.de@thermofisher.com Helpdesk E-Mail: service.lpg.germany.de@ThermoFisher.com

Anfragen aus Europa, mittlerem Osten und Afrika:

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

Postanschrift 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 Emai: linfo.labequipment@thermofisher.com

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

Enquiries from Asia Pacific: Phone +852-2711 3910 Fax +852-2711 3858 Emai: linfo.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

Anfragen von Europa:

Österreich Vertrieb:+43 1 801 40 0 Service:+43 1 801 40 0

Belgien Vertrieb:+32 53 73 4241 Service:+32 53 73 4241

Finnland/ Nordische/ Baltische Länder

Vertrieb:+358 9 329 100 Service:+358 9 329 100

Frankreich

Vertrieb:+33 2 2803 2180 Service:+33 825 800 119 Deutschland: Postanschrift Deutschland: Thermo Electron LED GmbH Robert-Bosch-Straße 1 D - 63505 Langenselbold

Telefon

Vertrieb Gebührenfrei 0800 1 536 376 oder +49 6184 90 6940 Service Gebührenfrei 0800 1 112110 oder +49 6184 90 6940 E-Mail info.labequipment.de@thermofisher.com

Italien

Vertrieb +39 02 95059 341 Service+39 02 95059 250

Niederlande

Vertrieb +31 76 579 5555 Service+31 76 579 5639

Russland/GUS

Vertrieb +7 812 703 4215 Service+7 812 703 4215

Spanien/Portugal

Vertrieb +34 93 223 0918 Service+34 93 223 0918

Schweiz Vertrieb +41 61 716 7755 Service+41 61 716 7755

England/ Irland Service+44 870 609 9203 Vertrieb+44 870 609 9203

[©] 2013 Thermo Fisher Scientific Inc. Alle Rechte vorbehalten. Alle Marken sind Eigentum von Thermo Fisher Scientific Inc. und ihrer Tochtergesellschaften. Spezifikationen, Bedingungen und Preise sind freibleibend. Nicht alle Produkte sind in allen Ländern verfügbar. Genauere Informationen sind auf Anfrage bei Ihrem lokalen Vertriebspartner erhältlich.

14 Kontaktdaten Thermo Scientific

6-fach-Gasblende für Cell Locker

Die 6-fach-Gasblende für Cell Locker wurde als segmentierte Innentür für das Cell Locker-System konzipiert.

Die 6-fach-Gasblende für Cell Locker unterteilt den vorhandenen Innenraum des CO_2 -Inkubators in sechs einzelne Nutzräume (Cell Locker). Zur Entnahme von Proben sind in der Frontpartie Glastüren (Zugangsöffnungen) eingelassen. Diese können separat geöffnet und geschlossen werden.



6-fach-Gasblende für CELL LOCKER



CELL LOCKER

Hinweis Geänderte Geräteleistung Bedingt durch die Ausführung der sechstürigen Gasblende ändern sich die Leistungsdaten des Inkubators (siehe <u>Technische Daten</u>).

Einlagebleche einschieben



Die Auflagebügel (1) in die unteren, mittleren und oberen **Vierkantlöcher** der Tragprofile einsetzen. Oben und in der Mitte die Lochblecheinlagen mit durchgängigen Führungsschienen einsetzen.



Abbildung: Lochblecheinlage mit durchgängigen Führungsschienen

Silikonstopfen nicht entfernen (1).

Bevor die Lochblecheinlage mit durchgängigen Führungsschienen platziert werden kann, müssen die Führungsschienen (2) eingeschoben werden.

Die Führungsschienen können ggf. entfernt werden.

Die Lochblecheinlage mit den durchgängigen Führungsschienen (1) und mittigen Öffnungen unten einsetzen.



Abbildung: Lochblecheinlage mit geteilten Führungsschienen für die untere Position

Diese Führungsschienen können nicht entfernt werden.



Abbildung: Einsetzen der Lochblecheinlage

Die Lochblecheinlagen so einsetzen, dass sie hinter der Schiene einrastet. Die Blechzunge an der Unterseite der Einlage muss hinter der Schiene sitzen, damit die Tür schließt.

Die 6-fach-Gasblende schließen.

A 6-fach-Gasblende für Cell Locker Einlagebleche einschieben

Betriebsanleitung

Cell Locker

50151663 Version März 2018



Vorwort

© 2018 Thermo Fisher Scientific Inc. Alle Rechte vorbehalten.

Thermo Fisher Scientific Inc. stellt seinen Kunden dieses Dokument nach Erwerb eines Produktes für den Betrieb des Gerätes zur Verfügung. Dieses Dokument ist urheberrechtlich geschützt. Die Vervielfältigung - auch auszugsweise – ist ohne schriftliche Zustimmung von Thermo Fisher Scientific Inc. verboten.

Änderungen an den Inhalten dieses Dokuments bleiben auch ohne Vorankündigung jederzeit vorbehalten.

Die in diesem Dokument enthaltenen Systemkonfigurationen und technischen Daten ersetzen etwaige frühere Angaben, die der Käufer erhalten hat.

Thermo Fisher Scientific Inc. erhebt keinen Anspruch auf die Vollständigkeit, Korrektheit und Fehlerfreiheit dieses Dokuments und haftet weder für hierin möglicherweise enthaltene Fehler oder Auslassungen noch für Folgeschäden, die sich aus der Verwendung dieses Dokuments ergeben, selbst wenn diese aufgrund der in diesem Dokument enthaltenen Angaben erfolgen sollten.

Dieses Dokument ist nicht Bestandteil eines Kaufvertrages zwischen Thermo Fisher Scientific Inc. und einem Käufer. Dieses Dokument hat keinerlei Änderungseinfluss auf die Allgemeinen Verkaufsbedingungen, vielmehr haben die Allgemeinen Verkaufsbedingungen bei voneinander abweichenden Angaben in den Dokumenten in jedem Fall Vorrang.

Für Übersetzungen in Fremdsprachen ist die deutsche Fassung dieser Betriebsanleitung verbindlich.

Warenzeichen

CELL LOCKER® ist ein eingetragenes Warenzeichen von Thermo Scientific. Thermo Scientific ist eine Marke der Thermo Fisher Scientific Inc.

Alle anderen in der Betriebsanleitung genannten Marken sind ausschließliches Eigentum der betreffenden Hersteller.

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

Die Thermo Electron LED GmbH ist eine Tochtergesellschaft von:

Thermo Fisher Scientific Inc. 168 Third Avenue Waltham, MA 02451 USA
Inhaltsverzeichnis

Kapitel 1 Allgemeines 2
1.1. Erklärung der Sicherheitshinweise und Bildzeichen 2
1.2. Sicherheitshinweise 3
1.3. Hinweise für den sicheren Betrieb 3
1.4. Gewährleistung 3
1.5. Verwendungszweck 4
1.5.1. Bestimmungsgemäßer Gebrauch 4
1.5.2. Bestimmungswidriger Gebrauch 4
Kapitel 2 Lieferumfang 5
Kapitel 3 Gerätebeschreibung
3.1. Beschreibung
3.2. Handhabung
3.3. Applikationsversionen 7
3.4. Erstverwendung
Kapitel 4 Reinigung und Desinfektion
4.1. Reinigung
4.2. Wisch- und Sprühdesinfektion 10
4.3. Autoklavierung
Kanitel 5 Service 11
5.1 Membran Wechsel
5.2. Dichtungswechsel Abdeckung
y 2 Prentango weekser ins deen ang i thirt i thirt i thirt i thirt i the
Kapitel 6 Ersatz- und Verschleißteile 14
Kapitel 7 Entsorgung 15
Kapitel 8 Technische Daten 16

1. Allgemeines

1.1 Erklärung der Sicherheitshinweise und Bildzeichen



Warnung

Bezeichnet eine Gefahrensituation, die zu Tod oder schwerer Verletzung führen könnte, wenn sie nicht vermieden wird!



Vorsicht

Bezeichnet eine Gefahrensituation, die zu leichter oder mäßiger Verletzung führen könnte, wenn sie nicht vermieden wird.

Achtung

Bezeichnet eine Situation, die zu Sachschäden führen könnte, wenn sie nicht vermieden wird.

	Sicherheitsrelevante Kapitel und Abschnitte innerhalb der Betriebsanleitung sind mit diesem Zeichen gekennzeichnet. Am Gerät angebracht verweist dieses Zeichen auf die besondere Beachtung der Betriebsanleitung, bzw. der Begleitpapiere.
T	Hinweis zur optimalen Nutzung des Gerätes
\bigtriangleup	Gefährliche Flüssigkeiten!
	Feuergefahr!
	Giftige Stoffe!
	Schutzhandschuhe tragen!
	Schutzbrille tragen!
	Mund- und Nasenschutz tragen!

1.2 Sicherheitshinweise

Der CELL LOCKER darf nur durch unterwiesenes Personal verwendet werden.

Für Personal, das mit dem CELL LOCKER arbeitet, sind auf der Grundlage dieser Betriebsanleitung, der gültigen Sicherheitsdatenblätter, der betrieblichen Hygienerichtlinien und der entsprechenden Technischen Regeln vom Betreiber schriftliche Verfahrensanweisungen zu erstellen, insbesondere:

- welche Dekontaminationsmaßnahmen für den CELL LOCKER und die verwendeten Hilfsmittel anzuwenden sind,
- welche Schutzmaßnahmen bei der Bearbeitung bestimmter Agenzien einzuhalten sind,
- welche Maßnahmen bei Unfällen zu ergreifen sind,
- zum Tragen von Schutzausrüstung, z. B beim Umgang mit mikrobiologischen und biologischen Proben.
- Trocknungsdauer 2 Stunden nach Außerbetriebnahme, z. B. wegen Transport.
- Arbeiten nach GLP Gute Laborpraxis.

1.3 Hinweise für den sicheren Betrieb

Wird der CELL LOCKER in ein Gerät eingebracht, ist dessen bestimmungsgemäßer Gebrauch und zulässige Gesamtlast zu beachten.

Die Dichtung der Abdeckung ist 1/2-jährlich auf Funktion und Schäden zu überprüfen.

Es dürfen keine Proben mit gesundheitsgefährlichen chemischen Substanzen verarbeitet werden, die durch Undichtigkeiten des CELL LOCKER in die Umgebungsluft freigesetzt werden oder auf Teile des Laborgerätes korrodierend, bzw. auf andere Art und Weise schädigend wirken können.

1.4 Gewährleistung

Thermo Fisher Scientific gewährleistet die Sicherheit und die Funktionstüchtigkeit des CELL LOCKER, für 1 Jahr, nur unter der Bedingung, dass:

- der CELL LOCKER ausschließlich für den bestimmungsgemäßen Gebrauch eingesetzt und gemäß den Angaben dieser Betriebsanleitung bedient und instand gehalten wird,
- keine baulichen Veränderungen am Gerät vorgenommen werden,
- nur originale und von Thermo Fisher Scientific zugelassene Ersatzteile oder Zubehöre verwendet werden (die Verwendung von Ersatzteilen von Fremdanbietern ohne Zulassung durch Thermo Fisher Scientific führt zum Erlöschen der Gewährleistungsansprüche),
- Inspektionen und Wartungsarbeiten entsprechend den vorgegebenen Zeitintervallen durchgeführt werden,
- nach allen Instandsetzungsarbeiten eine Funktionsprüfung durchgeführt wird.

Der Gewährleistungszeitraum beginnt mit der Auslieferung des CELL LOCKER an den Betreiber.

1.5 Verwendungszweck

1.5.1. Bestimmungsgemäßer Gebrauch

Der CELL LOCKER ist zur Verwendung in Laborgeräten vorgesehen, die unter anderem zur Aufbereitung und Kultivierung von Zell- und Gewebekulturen dienen.

Die maximale Autoklavier-Temperatur beträgt 121°C (250 °F).

1.5.2. Bestimmungswidriger Gebrauch

Keine Zell- oder Gewebekulturen im Gerät verwenden, die nicht den Vorschriften der Sicherheitsstufen L1, L2 und L3 entsprechen.

Es dürfen keine Gewebe, Stoffe oder Flüssigkeiten als Proben verwendet werden, die:

- leicht entflammbar oder explosiv sind,
- deren Dämpfe in Verbindung mit Luft brennbare oder explosive Gemische bilden,
- die Gifte freisetzen,
- Stäube bilden,
- exotherme Reaktionen verursachen,
- pyrotechnische Substanzen sind.

Außerdem dürfen keine Flüssigkeiten auf den Innenboden gegossen oder in einer Wanne in den Innenraum eingebracht werden.

Der Cell Locker darf nicht in Medizingeräten verwendet werden!

2. Lieferumfang

Beschreibung	Anzahl
Membrankammer	1
Satz Membranrahmen	1
Arbeitsplatte mit Abdeckung	1
Kleinteilesatz	1
Bedienungsanleitung	1
Zusammengefasste Sicherheitshinweise	1

Lieferkontrolle

Überprüfen Sie sofort nach Anlieferung des Gerätes:

- die Vollständigkeit der Lieferung,
- den Lieferzustand des Gerätes.

Ist die Lieferung unvollständig oder sind Transportschäden an Gerät und Verpackung feststellbar, informieren Sie umgehend den Spediteur, sowie den technischen Support von Thermo Fisher Scientific.

3. Gerätebeschreibung

3.1 Beschreibung

Der CELL LOCKER besteht aus einer Polycarbonatkammer mit zwei Membranfiltern, die für Wasserdampf und Gas durchlässig, aber für Mikroorganismen undurchlässig sind.



Abb. 1. Cell Locker

Position	Bezeichnung	
1	Arbeitsplatte	
2	Membranrahmen mit Filter	
3	Dichtung	
4	Verschluss	
5	Abdeckung	
6	Griff	

Produktionsjahr Produkti-

Polycarbonat Makrolon 2528

Abb. 2. Bodenmarkierungen

3.2 Erstverwendung

Vor dem ersten Gebrauch den CELL LOCKER reinigen und desinfizieren.



Hinweis

CELL LOCKER vor dem Einsetzen der Membranen reinigen.

3.3 Handhabung

Der CELL LOCKER kann mit einer gasdichten 6-türigen Gasblende mit oder ohne Abdeckung in den Inkubator gestellt werden.

Beim Einsetzen in ein Laborgerät darauf achten, dass der CELL LOCKER nicht verkantet.



3.4 Applikationsversionen

Je nach Kundenwunsch kann der CELL LOCKER mit einem Kleinteilesatz modifiziert werden, um die Abdeckung anzubringen oder zu entfernen.:



Abb. 3. Kleinteilesatz

Die Abdeckung von der Arbeitsplatte kann mit einem Inbusschlüssel entfernt werden, so dass durch die einzelnen Innentüren auf die Proben zugriffen werden kann.



mit installierter Abdeckung



ohne Abdeckung

Abb. 4. CELL LOCKER Optionen

4. Reinigung und Desinfektion

Vor Beginn der Arbeiten sind die folgenden Sicherheitshinweise zu beachten:



Warnung

Gesundheitsgefährdung!



Die Oberflächen des Nutzraums können kontaminiert sein. Der Kontakt mit kontaminierten Reinigungsflüssigkeiten kann Infektionen verursachen. Desinfektionsmittel können gesundheitsschädliche Stoffe enthalten.



Schutzbrille tragen!



Sicherheitshandschuhe tragen!



Mund- und Nasenschutz tragen!

Hinweise des Desinfektionsmittelherstellers und der Hygienefachkraft beachten!

4.1 Reinigung



Vorsicht

Unverträgliche Reinigungsmittel!

Teile des CELL LOCKER sind aus Kunststoff gefertigt.

Lösungsmittel können Kunststoffe anlösen.

Starke Säuren oder Laugen können eine Versprödung der Kunststoffe verursachen.



Vorsicht

Feuchtigkeitsempfindliche Bauteile!

Den Membranrahmen nicht mit Reinigungsmitteln besprühen. Beim Abwischen darauf achten, dass keine Feuchtigkeit eindringt.

Schmutzrückstände und Ablagerungen mit lauwarmem Wasser, das mit handelsüblichem Spülmittel versetzt ist, gründlich beseitigen.

Die Oberflächen und Arbeitsplatte mit einem sauberen Tuch und klarem Wasser abwischen.

Abschließend die Oberflächen mit einem sauberen Tuch trocken reiben.

4.2 Wisch- und Sprühdesinfektion



Vorsicht

Alkoholische Desinfektionsmittel!

Desinfektionsmittel, die mehr als 10 % Alkohol enthalten, können zusammen mit Luft leicht entflammbare und explosive Gasgemische bilden.

Bei Anwendung solcher Desinfektionsmittel offenes Feuer oder starke Hitzeeinwirkung während des gesamten Desinfektionsverfahrens vermeiden!

Solche Desinfektionsmittel nur in gut belüfteten Räumen anwenden.

Nach Einwirkung des Desinfektionsmittels die behandelten *CELL LOCKER* Geräteteile gut trocken reiben.

Sicherheitsregeln zur Vermeidung von Brand- und Explosionsgefahren durch alkoholische Desinfektionsmittel beachten.



Vorsicht

Chloridhaltige Reinigungsmittel!

Keine chloridhaltigen Desinfektionsmittel verwenden!

Zur Desinfektion nur Desinfektionsmittel benutzen, die sich unschädlich auf Edelstahl auswirken!



Warnung

Dekontaminations- oder Reinigungsmittel

Thermo Electron LED GmbH oder dessen Vertreter ist zu konsultieren, wenn irgendwelche Zweifel hinsichtlich der Verträglichkeit von Dekontaminatios- oder Reinigungsmitteln mit Teilen des Gerätes oder mit darin enthaltenen Stoffen bestehen.

- 1. Alle Proben aus dem Arbeitsbereich entfernen und an einem sicheren Ort aufbewahren.
- 2. Die Oberflächen mit Desinfektionsmittel abwischen. Desinfektionsmittel nicht auf die Membranfilter sprühen.
- 3. Warten, bis das Desinfektionsmittel die vom Hersteller angegebenen Werte erreicht hat.
- 4. Die Kammerflächen und die Arbeitsfläche mit lauwarmem Wasser abwischen, das mit Standardspülmitteln vermischt ist. Spülen Sie vollständig mit warmem Wasser ab.
- 5. Die gereinigten Flächen 3 bis 5 Mal mit autoklaviertem Wasser abspülen, um eventuelle Reinigungsmittelreste vollständig zu entfernen.
- 6. Anschließend die gereinigten Oberflächen, das Luftleitblech und die Einbauten mit einem weichen, sterilen Tuch abwischen.

4.3 Autoklavierung

Der CELL LOCKER kann bis zu 12 mal bei einer Temperatur von 121 °C autoklaviert werden.



Hinweis Beim Cell Locker vor dem Autoklavieren die Membranen entfernen!.

5. Service

Servicearbeiten sollten dokumentiert werden.

5.1 Membran Wechsel

Periodisch die Membranen überprüfen und spätestens nach einem Jahr die Membranen wechseln.



Warnung Gesundheitsgefährdung!



Schutzhandschuhe tragen!



Mund- und Nasenschutz tragen!

1. Arbeitsplatte bis zum Anschlag herausziehen und entnehmen.



2. Membrane von innen nach außen drücken, entnehmen.





3. Die Membranen in eine Plastiktüte stecken, autoklavieren und als Sondermüll entsorgen.





4. Die neuen Membranen auspacken.



- 5. Neue Membran ansetzen und im Rundlauf fest eindrücken.
- 6. Sichtprüfung von Innenseite, ob Dichtlippe korrekt sitzt, ggf. von außen nachdrücken.



7. Die zweite Membrane entsprechend Schritt 5 und 6 einsetzen.



Warnung Kontaminationsgefahr! Bei beschädigten oder falsch montierten Membranen kann der Anwender oder die Umgebung kontaminiert werden.!

8. Die Arbeitsplatte über Anschlag einsetzen und in Innenraum hineinschieben.



5.2 Dichtungswechsel Abdeckung

Periodisch, z.B. monatlich, die Dichtung der Abdeckung auf Intaktheit, keine Beschädigungen prüfen und falls erforderlich ersetzen.

1. Die Arbeitsplatte herausziehen und entnehmen.



2. Dichtung aus Nut entnehmen.



3. Neue Dichtung in Nut einsetzen und darauf achten, dass die Dichtungslippe nach außen zeigt.





WARNUNG Kontaminationsgefahr! Bei beschädigten oder falsch montierter Dichtung kann der Anwender oder die Umgebung kontaminiert werden.!

4. Die Arbeitsplatte in Innenraum hinein schieben.



6. Ersatz- und Verschleißteile

Beschreibung	Material-Nr.	
Membranrahmen (Satz mit 2)	50153148	
Dichtung Deckel	50151658	
Arbeitsplatte rostfreier Stahl	50151654	
Arbeitsplatte Kupfer	50154738	
Kleinteilesatz	50153097	
Deckel Membrankammer	50151656	

Es dürfen nur originale und von Thermo Fisher Scientific zugelassene Ersatzteile oder Zubehöre verwendet werden (die Verwendung von Ersatzteilen von Fremdanbietern ohne Zulassung durch Thermo Fisher Scientific führt zum Erlöschen der Gewährleistungsansprüche).

7. Entsorgung



Warnung Kontominationag

Kontaminationsgefahr!

Der Cell Locker konnte zur Be- und Verarbeitung von infektiösen Substanzen eingesetzt worden sein. Das Gerät oder Teile des Gerätes können deshalb kontaminiert sein. Alle Komponenten des Gerätes müssen vor der Entsorgung dekontaminiert werden!

Die Bauteile des Gerätes sind gründlich zu reinigen und anschließend abhängig vom Einsatzzweck entweder zu desinfizieren oder zu sterilisieren.

Dem Entsorgungsgut ist eine Unbedenklichkeitserklärung mit genauen Hinweisen über die Durchführung der Dekontaminationsmaßnahmen beizufügen.

Alle Komponenten, außer den Membran-Filtern, des CELL LOCKER können nach entsprechender Dekontaminierung der geregelten Entsorgung zugeführt werden. Die Membran-Filter müssen entsprechend den national gültigen Richtlinien für Sondermüll entsorgt werden.

Übersicht der verwendeten Materialien

Komponente	Material	
Cell Locker Gehäuse und Abdeckung	Polycarbonat, Makrolon 2858	
Arbeitsplatte rostfreier Stahl	1.4301+2B / Stahl 10088-2	
Arbeitsplatte Kupfer	Kupfer	
Membranrahmen	ABS	
Membran-Filter	PES Supor-200R Membrane 0,2 µm PALL	

8. Technische Daten

Bezeichnung	Einheit	Wert
Außenabmessungen (B x H x T)	mm	185 x 156 x 380
Innenraumvolumen	I	ca. 5.4
Gewicht	kg	1,8
Flächenlast Arbeitsplatte	kg	5
Temperaturbereich	°C / °F	4 - 55 / 39 - 131
Autoklavier-Zyklen		max. 12
Autoklavier-Temperatur	°C / °F	121 / 250



Find out more at thermofisher.com