Séries 8000DH

Incubateur à CO₂ **à jaquette d'air avec cycle de décontamination** Manuel d'utilisation et d'entretien 7013540 Rév. 10





Modèle	Capteur*	Tension**
3540	CT	115
3541	CT	230
3542	IR	115
3543	IR	230

*CT désigne un capteur à conductivité thermique.

IR désigne un capteur à infrarouge.

**Toutes les unités fonctionnent sur courant 50/60 Hz.



Important Lisez ce manuel d'instructions. Si les instructions de ce manuel ne sont pas lues, comprises et respectées, l'unité peut être endommagée, les opérateurs blessés et les performances réduites. ▲

Avertissement Tous les réglages et les tâches de maintenance doivent être effectués par du personnel d'entretien qualifié. ▲



- Utiliser ce produit uniquement de la manière décrite dans la documentation du produit et dans ce manuel. Avant de l'utiliser, vérifiez que ce produit est adapté à l'usage prévu.
- Ne pas modifier les composants du système, en particulier le contrôleur. Utilisez OEM équipement de remplacement exact ou des pièces. Avant l'utilisation, vérifiez que le produit n'a pas été modifié d'aucune façon.
- Débranchez l'appareil de toutes les sources d'alimentation avant le nettoyage, le dépannage, ou d'effectuer tout autre entretien sur le produit ou ses contrôles. Pour déconnecter l'alimentation à l'incubateur, débranchez le cordon d'alimentation à l'arrière de l'incubateur. Notez que l'activation du commutateur à clé sur le panneau de commande avant sur la position Off ne suffit pas de couper le courant.

Avertissement L'utilisateur est responsable de l'exécution des procédures de décontamination appropriées lorsque des matières dangereuses sont déversées sur ou à l'intérieur de l'incubateur.

Les informations figurant dans ce manuel sont uniquement fournies à titre indicatif. Son contenu et le produit qu'il décrit peuvent être modifiés sans préavis. Thermo Scientific ne fait aucune déclaration et n'accorde aucune garantie relatives à ce manuel. Thermo ne sera en aucun cas tenu responsable des dommages directs ou fortuits produits par ou liés à l'utilisation de ce manuel.

Avertissement Si l'incubateur n'est pas utilisé de la manière spécifiée dans ce manuel d'utilisation, la protection fournie par la conception de l'équipement pourrait être altérée.

© 2023 Thermo Fisher Scientific. Tous droits réservés.



Instructions importantes d'utilisation et de maintenance. Lisez attentivement le texte d'accompagnement.



Risques électriques potentiels. Seules des personnes qualifiées doivent effectuer les procédures associées à ce symbole.



Lors des tâches d'entretien ou de réparation, l'équipement doit être placé hors tension et verrouillé pour éviter toute blessure.



Avertissement de risque d'asphyxie. Des concentrations élevées de CO2 et de N2 peuvent déplacer l'oxygène et provoquer asphyxie!



Levage Avertissement de danger. L'incubateur pèse plus de 200 livres (91kgs). Prendre des mesures de sécurité adéquates lors du déplacement de ce dispositif.



Marquage de l'équipement électrique et électronique, qui s'applique à l'équipement électrique et électronique relevant de la Directive 2012/19/EU (DEEE) et à l'équipement mis sur le marché après le 13 août 2005.

Ce produit a l'obligation de satisfaire aux prescriptions de la directive européenne 2012/19/EU relative aux déchets d'équipements électriques (DEEE). Il est marqué du symbole DEEE. Thermo Scientific a signé un contrat avec une ou plusieurs entreprises de recyclage/élimination des déchets dans chaque État membre de l'UE et ce produit doit être mis au rebut ou recyclé par leur intermédiaire. Des informations supplémentaires sur la conformité de Thermo à cette directive, la liste des entreprises de recyclage pour votre pays ainsi que des informations sur les produits Thermo seront mises à votre disposition sur le site www.thermofisher.com.

- ✓ Utilisez toujours l'équipement de protection approprié (vêtements, gants, lunettes de protection, etc.)
- ✔ Dissipez toujours le froid ou la chaleur extrêmes et portez des vêtements protecteurs.
- ✓ Suivez toujours les bonnes pratiques d'hygiène.
- ✓ Chaque individual est responsable de sa propre sécurité.



Notre service clients peut vous aider pour les questions commerciales et vous fournir des devis et des informations sur les tarifs. Nous pouvons enregistrer votre commande et vous renseigner sur les modalités de livraison ou demander à votre revendeur le plus proche de vous contacter. Tous nos produits sont répertoriés sur notre site Web et vous trouverez toutes nos coordonnées sur la page d'accueil.

Notre service clients peut vous aider pour les questions techniques et vous renseigner sur la configuration, le fonctionnement et le dépannage de otre matériel. Nous pouvons vous envoyer des pièces de rechange ou détachées ou planifier les déplacements sur site. Nous pouvons également vous renseigner sur l'extension de garantie disponible pour vos produits Thermo Scientifi

Quel que soit le ou les produits Thermo Scientific que ous utilisiez ou dont vous ayez besoin, n'hésitez pas à nous contacter pour parler avec nous de vos applications. Si vous rencontrez un problème technique, nous essaierons ensemble d'identifier le p oblème et vous aiderons à le régler vous-même ...par téléphone, sans envoyer de technicien sur place.

En cas d'intervention plus complexe, nous dépêcherons l'un de nos techniciens qualifiés et formés ou une entreprise agréée pour procéder à la réparation sur place. Si l'intervention est couverte par la garantie, nous prendrons en charge les coûts de la réparation.

Remarques sur la garantie

Informations à connaître avant de demander une intervention dans le cadre de la garantie

- Localisez les numéros de modèle et de série. Une étiquette signalétique est apposée sur l'unité ellemême.
- Pour toute réparation ou maintenance de l'équipement, ou pour toute demande technique ou concernant une application spéciale, contactez votre distributeur local.

Réparations NON couvertes par la garantie

- Etalonnage des paramètres de commande. Les étalonnages nominaux sont effectués à l'usine : typiquement ±1 °C pour la température, ±1 % pour les gaz et ±5 % pour l'humidité. Notre service après-vente peut fournir des étalonnages précis sur votre site. Cette prestation est facturée. L'étalonnage effectué après une réparation dans le cadre de la garantie est couvert par la garantie.
- Tout dommage résultant d'une eau de mauvaise qualité, de substances chimiques ou de produits de nettoyage caustiques pour les matériaux de l'équipement.
- Demandes de dépannage résultant d'une installation ou d'instructions de fonctionnement incorrectes. Toutes les corrections apportées aux éléments suivants seront facturées :
 - 1) Connexion au réseau électrique
 - 2) Branchements des tubes
 - 3) Détendeurs de gaz
 - 4) Réservoirs de gaz
 - 5) Mise à l'horizontale de l'unité
 - 6) Ventilation de la pièce
 - 7) Fluctuations indésirables de la température ambiante
 - 8) Toute réparation externe à l'unité
- Dommage résultant d'un accident, d'une altération, d'un mauvais usage, d'un mauvais traitement, d'un incendie, d'une inondation, d'une catastrophe naturelle ou d'une installation incorrecte.
- Réparations de pièces du système résultant de modifications non autorisées de l'unité.
- Tous coûts de main-d'œuvre autres que ceux spécifiés pendant la période de garantie pièces et main-d'œuvre, pouvant inclure une garantie supplémentaire sur les capteurs de CO₂, les moteurs de ventilateurs, les jaquettes d'eau, etc.

Table des matières

Section 1	Installation et démarrage	1-1
	Utilisation du clavier	1-3
	Blocs d'affichage	1-4
	Installation de l'incubateur	1-4
	Superposition des incubateurs	1-5
	Nettovage préliminaire	1-6
	Installation des étagères	1-7
	Filtre du port d'accès et couvercle du capteur de $CO_2 \dots$	1-8
	Installation du filtre d'échantillonnage de l'air	1-8
	Installation du filtre HEPA	1-9
	Mise à niveau de l'unité	1-9
	Branchement électrique	1-9
	Remplissage du bac humidificateur	.1-10
	Branchement sur l'alimentation en CO ₂	.1-11
	Démarrage de l'incubateur	.1-12
Section 2	Étalonnage	2-1
	Température	2-1
	Mesure du CO ₂ par infrarouge	2-2
	Mesure par Infrarouge du CO ₂	2-3
	Humidité relative	2-4

Section 3	Configuration	3-1
	Activation/Désactivation de l'alarme sonore	3-1
	Nouveau filtre HEPA	3-1
	Réglage du rappel REPLACE HEPA	3-2
	Réglage du code d'accès	3-2
	Réglage de la limite d'alarme basse de temp. (alarme de suivi) .	3-3
	Activation des contacts d'alarme basse température	3-3
	Réglage de la limite d'alarme basse de CO ₂ (alarme de suivi)	3-4
	Réglage de la limite d'alarme haute de CO2 (alarme de suivi)	3-4
	Activation des contacts d'alarme de CO ₂	3-5
	Réglage du zéro pour les capteurs CT de CO2	3-5
	Réglage du nouvel intervalle de mesure pour les capteurs CT	
	de CO ₂	3-6
	Réglage d'une limite d'alarme basse de HR	3-6
	Activation des contacts d'alarme de HR	3-7
	Activation de l'affichage Temp/HR	3-7
Section 4	Alarmes	4-1
	Panne du contrôleur de température TMP CNTR FRR	4-2
	Alarmes de défauts des capteurs	4-7
	Maintenance préventive	4-3
Section 5	Entretien courant	5-1
	Nettoyage de l'extérieur de l'incubateur	5-2
	Nettoyage des portes en verre	5-3
	Nettoyage du bac humidificateur	5-3
	Maintenance du filtre HEPA	5-3
	Remplacement des fusibles	5-3
	Remplacement du filtre d'échantillonnage de l'air	5-4
	Electronique	5-4
	Cycle de décontamination	5-8

Section 6	Options installées en usine6-1
	Report d'alarmes
Section 7	Caractéristiques

Section 1 Installation et démarrage



Figure 1-1. Composants

- Orifice d'échantillonnage des gaz de la cuve Utilisé pour échantillonner le contenu en CO₂ de la cuve à l'aide d'un Testeur CO₂ FYRITE ou d'un instrument similaire.
- Commutateur d'alimentation principal Mise hors tension.
- Panneau de commande Clavier, blocs d'affichage et indicateurs (Figure 1-2).
- Pieds de mise à niveau Servent à mettre l'unité à l'horizontale.
- Bouton du cycle de décontamination Sert à lancer le cycle de décontamination.

Remarque Les incubateurs peuvent être empilés les uns sur les autres. Informations plus bas. ▲ (DI) - Désactive l'alarme sonore (Section 4).

Voyant d'alarme - Clignote pendant un état d'alarme.

Commutateur de MODE - Permet de sélectionner les modes Run (Fonctionnement), Setpoints (points de consigne), Calibration (Étalonnage) et Configuration.

Bloc d'affichage des messages - Affiche l'état du système.

Voyants de sélection du mode:

RUN : menu de fonctionnement

SET : menu des points de consigne

CAL : menu Étalonnage

CON : menu de configuration

Voyant HEAT - S'allume lorsque les résistances sont activées.

Voyant INJ - S'allume pendant l'injection du CO₂ dans la cuve.

- **Bloc d'affichage T** (°C) Programmable pour afficher la température. Voir la section Configuration.
- **Bloc d'affichage RH** (%) Affiche le pourcentage d'humidité. Voir la section Configuration.

Bloc d'affichage CO² (%) - Programmable pour afficher le pourcentage de CO₂. Voir la section Configuration.

Touches fléchées (Scroll for Parameters) - Incrémentent ou décrémentent les valeurs numériques, ou basculent entre les différents modes.

ENTER - Enregistre la valeur sélectionnée en mémoire.



Niveau de CO2 en %

Figure 1-2. Panneau de commande

Utilisation du clavier

L'incubateur Series 8000DH peut être configuré dans quatre modes de base : Run (Fonctionnement), Setpoints (Points de consigne), Calibration et Configuration.

- Run est le mode par défaut pour un fonctionnement normal.
- Set permet de saisir des points de consigne pour le fonctionnement de l'incubateur.
- Calibration sert à étalonner les différents paramètres du système.
- Configuration permet de définir diverses options pour personnaliser le fonctionnement.

↓			
MODE	MODE)→ MODE	→ MODE
RUN	SETPOINT	CALIBRATION	CONFIGURATION
Default Mode (Mode par défaut)	Temperature (Température)	Temp Offset (Écart de tem- pérature)	Audible
	Overtemp (Surtemp)	CO ₂ Cal ¹	New HEPA Timer (Nouvelle tem- porisation HEPA)
	CO ₂	IR Cal ²	Replace HEPA Reminder (Rappel remplacement HEPA)
		RH Cal (Cal HR)	Access Code (Code d'accès)
			Temp Lo Alarm (Alarme basse de temp.)
			Temp Relay (Relais temp.)
			CO2 Lo Alarm (Alarme basse CO2)
			CO2 Hi Alarm (Alarme haute CO2)
			CO2 Relay (Relais CO2)
			CO2 Z & S #'s (zéro et intervalle de mesure du CO2) *
			RH Lo Alarm (Alarme basse de HR)
			RH Relay (Relais HR)
			Display Temp (Afficher temp.)
			Display RH (Afficher HR)
¹ Unités CT		Menus des unités	de base
² Unités IR		Menus des unités a	vec options
*Unités CT uniquemen	t		

Tableau 1-1. Options de chaque mode

Utilisation du clavier (suite)

- Flèches SCROLL FOR PARAMETERS : fait accéder l'opérateur aux paramètres des modes SET, CAL et CON. La flèche droite affiche le paramètre suivant, la flèche gauche le paramètre précédent.
- Flèches vers le haut/bas : incrémentent/décrémentent la valeur du paramètre sélectionné dans les modes SET, CAL et CON ou passent d'un paramètre à un autre.
- ENTER : cette touche doit être enfoncée pour l'enregistrement en mémoire de toutes les valeurs modifiées.
- Touche (): appuyez pour désactiver l'alarme sonore. Voir la section 4 pour les intervalles de rappel des alarmes.

Blocs d'affichage Affichage des messages : affiche en permanence l'état du système (Mode). Affiche "CLASS 100" ou "SYSTEM OK" en fonctionnement normal, ou des messages d'alarme si le système détecte une condition d'alarme (voir Section 4, Alarmes). Le message "CLASS 100" est un mécanisme de temporisation indiquant que, dans des conditions normales de fonctionnement et lorsque le filtre HEPA est installé, l'air à l'intérieur de la cuve remplit les normes de propreté de la classe 100 pour les particules de 0,5 micron ou plus par mètre cube d'air. Thermo met à votre disposition des informations plus détaillées sur la classification de la qualité de l'air Classe 100.

Blocs d'affichage supérieur et inférieur : le bloc supérieur à 7 segments indique la température et l'humidité relative. Le bloc inférieur affiche le CO₂ et l'O₂.

Installation de l'incubateur

- 1. Maintenir un dégagement de six pouces minimum derrière l'unité pour branchement électrique et de gaz. En outre, un espace libre minimal de ventilation trois pouces est nécessaire de chaque côté.
- 2. Placez l'unité sur une surface solide et plane capable de supporter son poids de 118 kg plus le poids échantillon et équipement utilisateur total estimé.
- 3. Repérez l'incubateur loin des portes et fenêtres extérieures que les changements de la température extérieure et le contact avec la lumière solaire directe peuvent affecter les fonctions anti-condensation de l'unité.
- 4. Positionner l'ouverture de la porte loin de chauffage à air pulsé et conduits de refroidissement car ceux-ci peuvent transporter la poussière, la saleté et d'autres contaminants dans la chambre d'incubation, ainsi que affecter négativement la porte d'ouverture de temps de récupération.
- 5. Soulevez l'unité par les côtés de la base de l'armoire. Ne pas essayer de le soulever par l'avant et l'arrière. Cela met l'accent sur les charnières de la porte extérieure.

Superposition des incubateurs

Avertissement Installez les unités empilées contre un mur ou une structure similaire. ▲

Avertissement Lorsque les incubateurs sont empilés, ne laissez pas les deux portes extérieures ouvertes en même temps. ▲

Avertissement Si les unités ont fonctionné, éteignez-les toutes les deux et débranchez-les du secteur avant d'entamer toute intervention de maintenance. ▲



Deux pattes d'empilement (représentées à gauche) sont fournies dans le sachet de pièces livré avec chaque incubateur.

Figure 1-3. Pattes d'empilement

- 1. Retirez à le couvercle protégeant le cordon de la porte. Voir Figure 1-4. Retirez la fiche du connecteur.
- 2. Retirez les quatre vis fixant les charnières de la porte à l'unité. Retirez la porte et posez-la à l'écart.





Figure 1-4. Couvercle

Figure 1-5. Emplacements des bouchons

- 3. Dévissez les deux bouchons sur la face supérieure de l'incubateur du bas (Figure 1-5).
- 4. Dévissez et retirez les 4 pieds de mise à niveau de l'unité à placer en haut et soulevez-la pour l'installer sur l'unité du bas. Alignez tous les côtés.

Avertissement Cet incubateur pèse 118 kg. Munissez-vous de suffisamment de personnel pour soulever l'unité. Soulevez l'unité par les côtés à la base du boîtier afin d'éviter d'exercer des contraintes sur les charnières de la porte extérieure. ▲



Figure 1-6. Orifices des vis

Superposition des incubateurs (suite)

- 5. Insérez les pattes d'empilement dans les fentes à l'arrière des unités à empiler comme indiqué Figure 1-7.
- 6. Alignez les fentes des pattes sur les orifices de montage à l'arrière de l'incubateur du haut. Fixez les pattes à l'aide des vis et des rondelles fournies dans le sachet de pièces détachées. Voir Figure 1-8.



Figure 1-7. Patte dans la fente Figure 1-8. Pattes installées à l'arrière de l'unité

 Enfilez une vis 1/4 x 20 et sa rondelle fournies avec les pattes d'empilement, dans l'orifice situé derrière le couvercle. Ne serrez pas. Reportez-vous Figure 1-9.



8. Retirez le couvercle de la même zone de l'autre côté de l'unité supérieure.

Figure 1-9. Orifice de vis

- 9. Enfilez l'autre vis $1/4 \ge 20$ et sa rondelle dans cet orifice.
- 10. Serrez les vis des deux côtés.
- 11. Assemblez les charnières de porte sur l'unité. Fixez-les à l'aide des vis.
- 12. Enfilez le cordon de la porte dans le connecteur, comme précédemment. Fixez le couvercle.
- 13. Installez le couvercle de l'autre côté de l'unité.
- 14. Les incubateurs empilés sont prêts à fonctionner.

Nettoyage préliminaire

- 1. Retirez les revêtements en vinyle des montants des étagères, des panneaux latéraux et du conduit d'air, le cas échéant.
- 2. À l'aide d'un désinfectant de laboratoire approprié, tel que quaternary ammonium, nettoyez toutes les surfaces intérieures.

Attention Avant d'utiliser toute méthode de nettoyage ou de décontamination, sauf celles qui sont recommandées par le fabricant, les utilisateurs doivent vérifier avec le fabricant que la méthode proposée ne sera pas endommager l'équipement. Les déversements accidentels de matières dangereuses sur ou à l'intérieur de cet appareil sont la responsabilité de l'utilisateur. ▲

Installation des étagères

- 1. Installé inclus oeillets dans la bride arrière de chaque feuille conduit; 6 œillets par feuille. Voir la Figure 1-10.
- 2 Installez les grands panneaux métalliques latéraux en orientant les languettes vers le centre de la cuve, les fentes vers le haut. Chaque panneau peut être indifféremment placé à gauche ou à droite. Il suffit d'en retourner un pour l'adapter au côté opposé. Inclinez les panneaux latéraux en les plaçant dans la cuve pour que leurs parties supérieures s'enfilent dans le conduit d'air supérieur, puis guidez-les pour les placer en



Figure 1-10. Oeillets sur le dos de la feuille de conduit

position verticale. La Figure 1-11 représente le panneaux latéral orienté pour le côté droit de la cuve.



Figure 1-11. Montants des étagères et Figure 1-12. Montants et fentes panneaux latéraux

- 3. En vous reportant à la Figure 1-11, observez que les montants d'étagères gauche et droit ne présentent entre eux aucune différence.
- 4. Installez les montants d'étagères en plaçant la fente arrière du montant sur la languette arrière appropriée du panneau latéral. Tirez le montant vers l'avant et engagez sa fente avant sur la patte appropriée du panneau latéral. Reportez-vous Figure 1-12.
- 5. La figure 1-11 représente l'un des montants installé sur le panneau latéral droit.

Installation des étagères (suite)

6. Localiser la tige fournie, le printemps et embouts, 4 chacun. Monter le ressort à la tige par le positionnement de la fin du printemps au cours de la crête, à la tige. Ensuite, appuyez sur le ressort de l'autre côté. Il doit s'enclencher. Voir les figures 1-13 et 1-14.



Figure 1-14. Printemps Monter à Rod

7. Embout Glisser sur ressort (Ffigure 1-15).

Figure 1-13. Fin du Printemps

Remarque Lorsque l'installation des tiges, l'extrémité du ressort peut être installé sur chaque côté de la chambre. La tige supérieure avant est élevée dans la chambre et la tige arrière supérieur est inférieure à permettre l'accès à la porte d'accès de port de filtre. A





- 8. Installer une extrémité de la tige dans le trou approprié dans la feuille de conduit.
- 9. Comprimer le ressort (sous le capuchon d'extrémité) pour insérer

l'autre extrémité de la tige dans le trou dans la feuille de conduit sur le côté opposé.

10. Figure 1-16 montre les quatre tiges installées.



Figure 1-16. Four Rods Installed in Chamber

Filtre du port d'accès et couvercle du capteur de CO₂

- 1. Repérez l'ouverture dans le coin supérieur gauche de la cuve intérieure. Retirez la bande adhésive située sur la face extérieure du port d'accès.
- 2. Repérez le bouchon avec filtre dans le sachet fourni avec le matériel. Soulevez le couvercle situé sur le port d'accès métallique et installez l'ensemble bouchon avec filtre dans le port d'accès, à l'intérieur de la cuve. Voir Figure 1-17.
- Le sachet fourni avec le matériel contient également le couvercle du capteur de CO₂. Installez-le à l'aide des fixations 1/4 de tour. Reportez-vous Figure 1-17 pour l'emplacement.



Figure 1-17. Filtre et couvercle

Installation du filtre d'échantillonnage de l'air

- 1. Retirez le filtre de son sachet.
- 2. Installez le filtre d'échantillonnage de l'air sur la cannelure du tuyau situé derrière le conduit supérieur.
- 3. Insérez l'autre extrémité de l'ensemble sur la tubulure métallique du conduit supérieur. Reportez-vous Figure 1-18.



Figure 1-18. Installation du filtre d'échantillonnage de l'air

Installation du filtre HEPA

Mise en garde Manipulez le filtre avec précaution. Le matériau du filtre peut être endommagé s'il est manipulé sans précaution. Pour éviter d'endommager l'incubateur, n'utilisez pas l'unité sans que le filtre HEPA ne soit en place. ▲

- 1. Retirez le filtre de son emballage.
- 2. Retirez le revêtement plastique du filtre en veillant à ne pas toucher le matériau du filtre.
- 3. Installez le filtre comme indiqué Figure 1-13. Pour de plus amples informations sur la maintenance du filtre HEPA, reportez-vous Section 5.

Mise à niveau de l'unité

Vérifiez que l'unité est à l'horizontale en plaçant un niveau à bulle sur l'une des étagères. Tournez l'écrou hexagonal dans le sens antihoraire pour allonger le pied ou dans le sens horaire pour le raccourcir. Mettez l'unité à niveau sur l'axe d'avant en arrière et de gauche à droite.

Branchement électrique

Voir l'étiquette de série sur le côté de l'appareil pour les spécifications électriques ou se référer aux schémas électriques à la fin de ce manuel..

Mise en garde Les valeurs nominales de la plaque signalétique sont basées sur la charge de courant pendant le cycle de décontamination. Le courant de fonctionnement normal est nettement inférieur. Vérifiez que le circuit électrique peut supporter la charge de courant du cycle de décontamination. Branchez l'incubateur à un circuit dédié relié à la terre. Le connecteur du cordon d'alimentation constitue le dispositif de déconnexion du secteur pour l'incubateur. Placez l'unité de sorte qu'elle puisse être facilement débranchée. ▲

Branchement électrique (suite)

Branchez le cordon d'alimentation fourni sur la borne d'entrée d'alimentation à l'arrière du boîtier (Figure 1-19), puis sur un circuit électrique dédié et relié à la terre.



Figure 1-19. Arrière de l'incubateur

Remplir la casserole d'humidité avec de l'eau distillée stérile à l'intérieur de ¹/₂ "de la partie supérieure avec 0,8 gallons (3 litres) d'eau. Si l'eau ne peut pas être stérilisé à 2% solution d'ammonium quaternaire peut être utilisé comme un additif désinfectant contrôler les bactéries et d'autres contaminants organiques, le cas échéant. Placez la casserole directement dans le centre de la surface de l'incubateur pour assurer une humidité optimale et la réponse à la température.

Pour un fonctionnement optimal de l'incubateur, le type d'eau recommandée est stérile distillée type 2 l'eau avec un pH compris entre 7 et 9, et une résistivité entre 50k Ω - 1M Ω cm (conductivité de 1-20 μ S / cm). Osmose inverse (RO), désionisée (DI), et ultrapure Type 1 'eau peut corroder inoxydable acier, cuivre, verre, et d'autres composants incubateur au fil du temps en raison l'eau de haute pureté avec un contenu ionique basse et haute résistivité sera activement tirer des ions à partir de ces matériaux. Si distillée de type 2 de l'eau stérile est pas facilement disponibles, ces autres types d'eau peuvent être utilisés en ajoutant le quantité nécessaire d'une solution stérile de bicarbonate de sodium faible pour élever le pH entre 7 et 9 et abaisser la résistivité entre 50k Ω - 1M Ω -cm (conductivité de 1-20 μ S / cm).

Les systèmes de distillation, ainsi que certains types de pureté de l'eau par osmose inverse systèmes, peuvent produire de l'eau dans la gamme de qualité spécifiée. L'eau du robinet est pas recommandé car il peut contenir du chlore, qui peut détériorer la acier inoxydable. L'eau du robinet peut également avoir une teneur élevée en minéraux, qui produirait une accumulation de tartre dans la casserole. Pureté élevée ou ultra pur l'eau ne soit pas recommandé, car il est un solvant extrêmement agressif et détériorer l'acier inoxydable.

Remplissage du bac humidificateur

Remplissage du bac humidificateur (suite)

Branchement sur l'alimentation en CO₂



L'eau de haute pureté a une résistance supérieure à 1M à 18M Ohm. Même l'eau pureté élevée peut contenir des bactéries et des contaminants organiques. L'eau doit toujours être stérilisée ou traitée avec un décontaminant sans danger pour l'inox et pour le produit avant d'être introduite dans le bac humidificateur.

Vérifiez le niveau et changez l'eau fréquemment pour éviter toute contamination. Évitez toute variation significative de niveau. Les dessèchements peuvent affecter de façon négative le niveau d'humidité et l'étalonnage de CO₂ des capteurs CT.

Mise en garde L'eau distillée ou désionisée utilisée dans la casserole d'humidité doit être dans une plage de résistance à la qualité de l'eau de 50K à 1M Ohm / cm pour protéger et prolonger la durée de vie de l'acier inoxydable. Utilisation de l'eau en dehors de la spécifiée gamme va diminuer la durée de fonctionnement de l'appareil et peut annuler la garantie. ▲

Avertissement De hautes concentrations de CO_2 peuvent provoquer l'asphyxie ! Les normes OSHA stipulent que l'exposition du personnel au dioxyde de carbone pendant une vacation de huit heures dans une semaine de travail de 40 heures ne doit pas excéder la moyenne pondérée sur huit heures de 5000 ppm (0,5 % CO₂). La limite d'exposition à court terme pour 15 minutes ou moins est de 30 000 ppm (3 % CO₂). Des moniteurs de dioxyde de carbone sont recommandés dans les zones confinées où les concentrations de CO₂ peuvent s'accumuler.

Avertissement Cet incubateur est conçu pour fonctionner avec du CO₂ uniquement. Tout branchement d'un gaz inflammable ou toxique peut produire des situations dangereuses.

Aucun gaz autre que le CO₂ ne doit être branché sur cet équipement. Les bouteilles de CO₂ comportent une étiquette UN1013 et sont équipées d'une valve de sortie CGA 320. Vérifiez que la bouteille de gaz comporte les libellés d'identification appropriés. La source de CO₂ branchée sur l'incubateur doit être de niveau industriel et pure à 99,5 %. N'utilisez pas de bouteilles de CO₂ équipées de siphon. Un siphon est utilisé pour extraire du CO₂ liquide de la bouteille, ce qui peut endommager le régulateur de pression. Assurezvous auprès de votre fournisseur de gaz que votre bouteille de CO₂ ne contient pas de siphon. Les bouteilles de gaz doivent être fixées au mur ou à un autre objet fixe pour les empêcher de basculer.

Un régulateur de pression du CO₂ à deux étages doit être installé sur la valve de sortie de la bouteille de gaz. La pression d'entrée de l'incubateur doit être maintenue à 15 psig (103,4 kPa) pour un bon fonctionnement du système de contrôle du CO₂. Un régulateur de pression du CO₂ à un étage ne maintient pas 15 psig (103,4 kPa) dans l'incubateur lorsque la pression à l'intérieur de la bouteille de CO₂ diminue ; c'est pourquoi un régulateur à deux étages est recommandé.

Branchement sur l'alimentation en CO₂ (suite)

Avertissement Si un CO₂ plus pur est souhaité à l'intérieur de l'incubateur (pureté supérieure à 99,5 %), le régulateur de pression doit être construit avec un diaphragme en inox, et la pureté du CO₂ doit être spécifiée par le fournisseur du gaz. Suivez les instructions du fabricant pour assurer une installation appropriée et sûre du régulateur de pression sur la bouteille de gaz.

Assurez-vous auprès du responsable de la sécurité de votre établissement que l'équipement est installé conformément aux codes et réglementations applicables localement.

L'alimentation en CO_2 doit être de qualité industrielle, pure à 99,5 %, et ne pas contenir de siphons. Installez un régulateur de pression à deux étages sur la sortie de la bouteille. Le manomètre de haute pression doit proposer une gamme de 0 à 2000 psig. Le manomètre de basse pression, à l'entrée de l'incubateur, doit proposer une gamme de 0 à 30 psig. La pression à l'entrée de l'incubateur doit être maintenue à 15 psig (103,4 kPa).

L'incubateur comporte des raccords dentelés à l'arrière du boîtier destinés au branchement de l'alimentation en gaz. Reportez-vous Figure 1-14. Le raccord est libellé "CO₂ Inlet #1 Tank". Vérifiez que les branchements sont fixés par des colliers. Vérifiez l'absence de fuites sur chaque raccord.

Démarrage de l'incubateur

L'incubateur étant correctement installé et sous tension, le bac humidificateur rempli et l'unité branchée sur l'alimentation en gaz, les points de consigne du système peuvent être saisis. Les points de consigne suivants peuvent être saisis en mode SET : Température, Surtempérature et CO₂. Pour entrer en mode SET, appuyez sur la touche MODE jusqu'à ce que le voyant SET s'allume. Appuyez sur les touches droite et/ou gauche jusqu'à ce que le paramètre approprié s'affiche sur le bloc d'affichage. Voir le Schéma 1-1 pour plus de détails.

Réglage de la température de fonctionnement

Tous les incubateurs series 8000DH ont une gamme de température de fonctionnement de 10 °C à 50 °C, selon la température ambiante. L'incubateur est expédié de l'usine avec un point de consigne de température défini à 10 °C. À ce réglage, toutes les résistances sont désactivées.

Réglage de la température de fonctionnement (suite)

Pour modifier le point de consigne de la température de fonctionnement :

- 1. Appuyez sur la touche MODE jusqu'à ce que le voyant SET s'allume.
- 2. Appuyez sur la flèche droite jusqu'à ce que "TEMP XX.XC" s'affiche sur le bloc d'affichage.
- 3. Appuyez sur la flèche vers le haut/bas jusqu'à ce que le point de consigne de température souhaité s'affiche.
- 4. Appuyez sur ENTER pour enregistrer le point de consigne.
- 5. Appuyez sur la touche MODE jusqu'à ce que le voyant RUN s'allume pour entrer en mode RUN, ou sur les flèches droite/gauche pour passer au paramètre suivant/précédent.

Réglage du point de consigne de surtempérature

Mise en garde Le système indépendant de surtempérature est conçu comme une sécurité pour protéger uniquement l'incubateur. Il n'est pas destiné à protéger les cultures de cellules ou un équipement stocké à l'intérieur de l'incubateur ni à limiter leur température maximale en cas de température excessive. ▲

Les incubateurs series 8000DH sont équipés d'un système de surveillance secondaire de la température destiné à contrôler la température de l'air à l'intérieur de l'appareil. Ce système est conçu comme un dispositif de sécurité désactivant toutes les résistances en cas de panne de contrôle de la température. Le contrôle de la température dans l'incubateur est réglé sur ±1° du point de consigne de surtempérature.

La surtempérature est réglée en usine (par défaut) à 40 °C. Néanmoins, elle peut être réglée jusqu'à 55 °C par paliers de 0,5 °C.

Si le point de consigne de la température ambiante de l'incubateur est réglé sur une valeur supérieure à de celui de surtempérature, ce dernier passe automatiquement à 1 °C au-dessus du point de consigne de la température. Il est recommandé de maintenir le point de surtempérature à 1 °C au-dessus de celui de la température de fonctionnement.

Réglage du point de consigne de surtempérature (suite)

Pour définir le point de consigne de surtempérature :

- 1. Appuyez sur la touche MODE jusqu'à ce que le voyant SET s'allume.
- 2. Appuyez sur la flèche droite jusqu'à ce que "OTEMP XX.XC" apparaisse sur le bloc d'affichage.
- 3. Appuyez sur la flèche vers le haut ou vers le bas jusqu'à ce que la valeur souhaitée s'affiche.
- 4. Appuyez sur ENTER pour enregistrer ce paramètre.
- 5. Appuyez sur la touche MODE jusqu'à ce que le voyant RUN s'allume, ou sur la flèche droite ou gauche pour passer au paramètre suivant ou précédent.

Réglage du point de consigne de CO2
Toutes les cellules CT de CO2 sont calibrées à l'usine à 37 °C, avec une humidité élevée et 10 % de CO2. Pour cette raison, si un point de consigne de température de 37 °C a été saisi, si le bac humidificateur a été rempli et si le contrôle du CO2 est réglé entre 0 et 10 % avec un capteur CT de CO2, le point de consigne du CO2 peut être immédiatement saisi. Sinon, il est important de laisser l'unité se stabiliser 12 heures au point de consigne de la température avant de saisir le point de consigne du CO2.

Sur les deux modèles d'incubateur, la gamme des points de consigne du CO_2 s'étend de 0,0 % à 20,0 %. L'incubateur est expédié de l'usine avec un point de consigne du CO_2 à 0,0 %. Avec cette valeur, toutes les commandes et alarmes de CO_2 sont désactivées. Pour modifier le point de consigne du CO_2 :

- 1. Appuyez sur la touche MODE jusqu'à ce que le voyant SET s'allume.
- 2. Appuyez sur la flèche de droite jusqu'à ce que "CO2 XX.X%" apparaisse sur le bloc d'affichage des messages.
- 3. Appuyez sur la flèche vers le haut/bas jusqu'à ce que le point de consigne de CO₂ souhaité s'affiche.
- 4. Appuyez sur ENTER pour enregistrer le point de consigne.
- 5. Appuyez sur la touche MODE jusqu'à ce que le voyant RUN s'allume si vous souhaitez passer en mode RUN ou sur la flèche vers la droite/gauche pour passer au paramètre suivant/précédent.





Section 2 Étalonnage

Lorsque l'unité est stabilisée, plusieurs systèmes doivent être étalonnés. En mode Calibration, la température de l'air ainsi que les niveaux de CO₂ et de HR peuvent être étalonnés avec des instruments de référence. Pour accéder au mode Calibration, appuyez sur la touche MODE jusqu'à ce que le voyant CAL s'allume. Appuyez sur la flèche vers la droite ou la gauche jusqu'à ce que le paramètre approprié apparaisse sur le bloc d'affichage des messages. Reportez-vous au Schéma 2-1 à la fin de cette section pour plus de détails.

La fréquence des étalonnages dépend de l'utilisation, des conditions ambiantes et de la précision requise. Une bonne pratique de laboratoire consiste en au moins un étalonnage annuel. Sur les nouvelles installations, tous les paramètres doivent être vérifiés après la période de stabilisation.

Avant l'étalonnage, l'utilisateur doit être conscient des fonctions suivantes du système. Lorsque l'unité est en mode Calibration, toutes les fonctions de commande du système sont à l'arrêt pour que l'unité reste stable. Les valeurs du système étalonné apparaissent sur le bloc d'affichage des messages. Si aucune touche n'est enfoncée pendant environ cinq minutes en mode Calibration, le système revient en mode RUN et les fonctions sont réactivées.

Mise en garde Avant tout étalonnage ou réglage de l'unité, il est impératif d'étalonner correctement tous les instruments de référence. ▲

Température Avant l'étalonnage, laissez la température de l'incubateur se stabiliser. Placez l'instrument étalonné au centre de la cuve. L'instrument doit se trouver dans le courant d'air et non contre l'étagère.

Périodes de stabilisation de la température

Démarrage - Laissez la température se stabiliser 12 heures dans l'incubateur avant de continuer.

En cours de fonctionnement - Laissez la température se stabiliser au moins 2 heures après l'affichage du point de consigne avant de continuer.

Température (suite) 1. Appuyez sur la touche MODE jusqu'à ce que le voyant CAL s'allume.

- 2. Appuyez sur la flèche droite jusqu'à ce que "TEMPCAL XX.X" apparaisse sur le bloc d'affichage des messages.
- 3. Appuyez sur la flèche vers le haut/bas jusqu'à ce que la valeur affichée corresponde à la lecture de l'instrument étalonné.
- 4. Appuyez sur ENTER pour enregistrer l'étalonnage.
- 5. Appuyez sur la touche MODE pour revenir en mode RUN ou sur la flèche vers la droite/gauche pour passer au paramètre suivant/précédent.

Mesure du CO₂ par conductivité thermique

Les modèles 3540 et 3541 comportent un capteur de CO₂ à conductivité thermique (CT). La conductivité thermique de l'atmosphère de l'incubateur n'est pas uniquement affectée par la quantité de CO₂ présent mais également par la température de l'air et la vapeur d'eau présente. Lors de la surveillance des effets du CO₂, la température de l'air et l'humidité absolue doivent donc être maintenues constantes pour que toute variation de la conductivité thermique reflète une variation de la concentration de CO₂.

Tout changement de température ou passage du niveau d'humidité élevée au niveau d'humidité ambiant de la pièce nécessite un réétalonnage de la commande de CO₂.

Certains capteurs CT de CO₂ connaissent une période de vieillissement, particulièrement sur les nouvelles installations. Dans ce cas, l'étalonnage doit être effectué toutes les semaines et le système ajusté en conséquence. Lorsque le capteur se stabilise, les vérifications peuvent devenir moins fréquentes.

Périodes de stabilisation du capteur CT de CO2

Démarrage - Le capteur de CO₂ a été calibré à l'usine sur 37 °C et une humidité élevée. Laissez la température ainsi que les niveaux d'humidité et de CO₂ se stabiliser au moins 12 heures dans la cuve avant de vérifier la concentration de CO₂ à l'aide d'un instrument indépendant.

En cours de fonctionnement - Vérifiez que les portes de la chambre sont fermées. Laissez l'atmosphère de la cuve se stabiliser au moins 2 heures après que les valeurs affichées de température et de CO₂ aient atteint leurs points de consigne.

Mesure du CO₂ par conductivité thermique (suite)

- 1. Veillez à ce que les périodes de stabilisation indiquées ci-dessus soient respectées.
- 2. Échantillonnez l'atmosphère de la cuve par le port d'échantillonnage à l'aide d'un instrument indépendant. Échantillonnez l'atmosphère au moins 3 fois afin de vous assurer de la précision de l'instrument.
- 3. Appuyez sur la touche MODE jusqu'à ce que le voyant CAL s'allume.
- 4. Appuyez sur la flèche de droite jusqu'à ce que "CO2 CAL XX.X" apparaisse sur le bloc d'affichage des messages.
- 5. Appuyez sur les flèches vers le haut/bas pour faire correspondre la valeur affichée à celle de l'instrument indépendant.
- 6. Appuyez sur ENTER pour enregistrer l'étalonnage.
- 7. Appuyez sur la touche MODE pour revenir en mode RUN ou sur les flèches droite ou gauche pour passer au paramètre suivant ou précédent.

Mesure du CO₂ par infrarouge

Les modèles 3542 et 3543 sont équipés d'un capteur de CO₂ à infrarouge (IR). Les capteurs de CO₂ à infrarouge ne sont pas affectés par la température et l'humidité de l'atmosphère de la chambre. Néanmoins, le détecteur de lumière du capteur est affecté par de grandes variations de température. Pour cette raison, une modification des points de consigne de la température peut nécessiter un réétalonnage du CO₂. Laissez à la température de la cuve le temps de se stabiliser avant de vérifier les concentrations de CO₂ à l'aide d'un instrument indépendant, particulièrement au démarrage.

Temps de stabilisation du capteur IR de CO2

Démarrage - Laissez la température et le CO₂ du boîtier se stabiliser au moins 12 heures avant de continuer.

En cours de fonctionnement - Laissez le CO₂ se stabiliser au moins 2 heures au point de consigne avant de continuer.

Mesure du CO2 par infrarouge (suite)	 Mesurez la concentration de CO2 dans la cuve par le port d'échantillonnage du gaz à l'aide d'un Fyrite ou d'un autre instrument indépendant. Plusieurs lectures sont nécessaires pour assurer la précision.
	2. Appuyez sur la touche MODE jusqu'à ce que le voyant CAL s'allume.
	3. Appuyez sur la flèche droite jusqu'à ce que "IR CAL XX.X" apparaisse sur le bloc d'affichage des messages.
	4. Appuyez sur la flèche vers le haut/bas jusqu'à ce que la valeur affichée corresponde à la lecture de l'instrument indépendant.
	5. Appuyez sur ENTER pour enregistrer l'étalonnage.
	6. Appuyez sur la touche MODE pour revenir en mode RUN.
Humidité relative	Les incubateurs modèles 3540 et 3541 sont équipés en option d'un capteur d'humidité relative à lecture directe. Il s'agit uniquement d'une lecture du niveau d'humidité relative de la cuve. Ils ne permettent pas de commander l'humidité relative à l'intérieur de l'incubateur.
	Temps de stabilisation de l'humidité relative
	Démarrage - Laissez l'humidité relative et la température de la cuve se stabiliser 12 heures avant de continuer.
	En cours de fonctionnement - Laissez l'humidité relative se stabiliser 2 heures après que la température affichée ait atteint le point de consigne de l'humidité relative avant de continuer.
	 Placez un instrument indépendant précis au centre de la cuve. Laissez l'humidité relative se stabiliser au moins 30 minutes.
	2. Appuyez sur la touche MODE jusqu'à ce que le voyant CAL s'allume.
	3. Appuyez sur la flèche droite jusqu'à ce que "RH XX%" apparaisse sur le bloc d'affichage des messages.
	 Appuyez sur la flèche vers le haut/bas jusqu'à ce que la valeur affichée corresponde à la lecture de l'instrument indépendant.
	5. Appuyez sur ENTER pour enregistrer l'étalonnage.
	6. Appuyez sur la touche MODE pour revenir en mode RUN.

Humidité relative (suite)

Si vous ne disposez d'aucun dispositif fiable de mesure de l'HR, l'écran peut être étalonné à un niveau standard:

- 1. Respectez les périodes de stabilisation de l'HR indiquées ci-dessus.
- 2. Avec un bac humidificateur plein et une température stable, l'humidité relative de la cuve est de 95 %.
- 3. En suivant les étapes 3 à 5 du réglage du capteur d'humidité relative, réglez l'affichage sur 95 %.

Cette méthode d'étalonnage présente normalement une précision à 5 % ou moins.

Tableau 2-1. Mode Calibration



Section 3 Configuration

Plusieurs fonctions disponibles en mode Configuration permettent de personnaliser la configuration de l'incubateur. Ces fonctions sont répertoriées et décrites ci-dessous. Toutes les fonctions peuvent ne pas être nécessaires dans toutes les applications, mais elles sont disponibles en cas de besoin. Pour entrer en mode Configuration, appuyez sur la touche MODE jusqu'à ce que le voyant CON s'allume. Appuyez sur les flèches droite et/ou gauche jusqu'à ce que le paramètre approprié s'affiche à l'écran des messages. Voir le Schéma 3-1 pour plus de détails.

Activation/ Désactivation de l'alarme sonore

L'alarme sonore peut être activée ou désactivée. Elle est activée par défaut en usine.

- 1. Appuyez sur la touche MODE jusqu'à ce que le voyant CON s'allume.
- 2. Appuyez sur la flèche droite jusqu'à ce que "AUDIBLE XXX" apparaisse sur le bloc d'affichage des messages.
- 3. Appuyez sur la flèche vers le haut/bas pour activer ou désactiver l'alarme sonore.
- 4. Appuyez sur ENTER pour enregistrer ce paramètre
- 5. Appuyez sur la touche MODE pour revenir en mode RUN ou sur la flèche vers la droite/gauche pour passer au paramètre suivant/précédent.

Nouveau filtre HEPA Lorsque le rappel "REPLACE HEPA" s'affiche et que l'alarme visuelle clignote, la durée spécifiée s'est écoulée et le filtre HEPA doit être remplacé. Pour effacer le message et réinitialiser la temporisation après avoir remplacé le filtre HEPA par un filtre neuf, procédez comme suit.

- 1. Appuyez sur la touche MODE jusqu'à ce que le voyant CON s'allume.
- 2. Appuyez sur la flèche droite jusqu'à ce que "NEW HEPA" s'affiche.
- 3. Appuyez sur ENTER pour relancer la temporisation et effacer le message d'alarme REPLACE HEPA.
- 4. Appuyez sur la touche MODE pour revenir en mode RUN.

Réglage du rappel REPLACE HEPA

Une temporisation de rappel pour le remplacement du filtre HEPA peut être définie sur une durée spécifique comprise entre 1 et 12 mois de fonctionnement réel de l'unité. Le temps pendant lequel l'unité est hors tension n'est pas pris en compte dans la temporisation. La durée par défaut est de 6 mois. Lorsque la durée fixée est écoulée, le message "REPLACE HEPA" s'affiche et l'alarme visuelle clignote. Pour configurer le rappel, procédez comme suit.

- 1. Appuyez sur la touche MODE jusqu'à ce que le voyant CON s'allume.
- 2. Appuyez sur la flèche droite jusqu'à ce que "REPL HEPA XX" s'affiche.
- 3. Appuyez sur la flèche vers le haut/bas pour sélectionner le nombre de mois souhaité.
- 4. Appuyez sur ENTER pour enregistrer ce nombre.
- 5. Appuyez sur la touche MODE pour revenir en mode RUN ou sur la flèche vers la droite/gauche pour passer au paramètre suivant/précédent.

Remarque Lorsque le rappel a été défini, vérifiez le temps restant en passant en mode Configuration, puis en appuyant sur la flèche droite jusqu'à ce que "NEW HEPA XXX" s'affiche. XXX est le nombre de jours restant avant la fin de la temporisation. Si, par exemple, vous avez choisi 12 mois dans le message "REPL HEPA XX", la valeur affichée dans le message NEW HEPA sera 365 jours. ▲

Réglage du code d'accès

Un code d'accès à 3 chiffres peut être saisi pour éviter que du personnel non autorisé ne modifie les points de consigne, l'étalonnage ou la configuration. Une valeur de 000 désactive le code d'accès. La valeur par défaut définie en usine est 000.

- 1. Appuyez sur la touche MODE jusqu'à ce que le voyant CON s'allume.
- 2. Appuyez sur la flèche droite jusqu'à ce que "ACC CODE XXX" apparaisse sur le bloc d'affichage des messages.
- 3. Appuyez sur la touche haut/bas pour modifier le code d'accès.
- 4. Appuyez sur ENTER pour enregistrer le code d'accès.
- 5. Appuyez sur la touche MODE pour revenir en mode RUN ou sur la flèche vers la droite/gauche pour passer au paramètre suivant/précédent.

Réglage de la limite d'alarme basse de temp. (alarme de suivi)

La limite d'alarme basse de température est la déviation par rapport au point de consigne de la température qui provoque une alarme basse de température. Cette alarme basse peut être définie entre 0,5 °C et 5 °C sous le point de consigne. Le réglage usine est de 1 °C sous le point de consigne. Un signe moins (-) sur le bloc d'affichage indique que la valeur de l'alarme est en dessous du point de consigne.

- 1. Appuyez sur la touche MODE jusqu'à ce que le voyant CON s'allume.
- 2. Appuyez sur la flèche droite jusqu'à ce que "TEMP LO -X.X" apparaisse sur le bloc d'affichage des messages.
- 3. Appuyez sur la touche haut/bas pour modifier la limite d'alarme basse de température.
- 4. Appuyez sur ENTER pour enregistrer la limite d'alarme basse de température.
- 5. Appuyez sur la touche MODE pour revenir en mode RUN ou sur la flèche vers la droite/gauche pour passer au paramètre suivant/précédent.

Activation des contacts d'alarme basse température

L'alarme basse température peut être programmée pour déclencher les contacts d'alarme à distance. C'est le cas si le paramètre est réglé sur ON. Un réglage sur OFF empêche l'alarme basse température de déclencher les contacts. Le réglage usine est ON.

- 1. Appuyez sur la touche MODE jusqu'à ce que le voyant CON s'allume.
- 2. Appuyez sur la flèche droite jusqu'à ce que "TMP RLY XXX" s'affiche.
- 3. Appuyez sur la flèche vers le haut/bas pour activer (ON) ou désactiver (OFF) ce déclenchement.
- 4. Appuyez sur ENTER pour enregistrer ce paramètre
- 5. Appuyez sur la touche MODE pour revenir en mode RUN ou sur la flèche vers la droite/gauche pour passer au paramètre suivant/précédent.

Réglage de la limite d'alarme basse de CO₂ (alarme de suivi)

La limite d'alarme basse de CO₂ est la déviation par rapport au point de consigne du CO₂ qui provoque une alarme basse de CO₂. Sa valeur peut être définie entre 0,5 % de CO₂ et 5,0 % de CO₂ en dessous du point de consigne. Le réglage usine est de 1,0 % de CO₂ en dessous du point de consigne. Un signe moins (-) sur le bloc d'affichage indique que la valeur de l'alarme est en dessous du point de consigne.

- 1. Appuyez sur la touche MODE jusqu'à ce que le voyant CON s'allume.
- 2. Appuyez sur la flèche droite jusqu'à ce que "CO2 LO -X.X" apparaisse sur le bloc d'affichage des messages.
- 3. Appuyez sur la touche haut/bas pour modifier la limite d'alarme basse de CO₂.
- 4. Appuyez sur ENTER pour enregistrer la limite d'alarme basse de CO2.
- 5. Appuyez sur la touche MODE pour revenir en mode RUN ou sur la flèche vers la droite/gauche pour passer au paramètre suivant/précédent.

Réglage de la limite d'alarme haute de CO₂ (alarme de suivi)

La limite d'alarme haute de CO₂ est la déviation par rapport au point de consigne du CO₂ qui déclenche une alarme haute de CO₂. Sa valeur peut être définie entre 0,5 % et 5,0 % au-dessus du point de consigne du CO₂. Le réglage usine est de 1,0 % de CO₂ au dessus du point de consigne.

- 1. Appuyez sur la touche MODE jusqu'à ce que le voyant CON s'allume.
- 2. Appuyez sur la flèche droite jusqu'à ce que "CO2 HI X.X" apparaisse sur le bloc d'affichage des messages.
- 3. Appuyez sur la touche haut/bas pour modifier la limite d'alarme haute de CO₂.
- 4. Appuyez sur ENTER pour enregistrer la limite d'alarme haute de CO2.
- 5. Appuyez sur la touche MODE pour revenir en mode RUN ou sur la flèche vers la droite/gauche pour passer au paramètre suivant/précédent.

Activation des contacts d'alarme de CO₂

Les alarmes haute et basse de CO₂ peuvent être programmées pour déclencher les contacts d'alarme à distance. Un réglage sur ON active cette fonction, un réglage sur OFF empêche les alarmes de CO₂ de déclencher les contacts. Le réglage usine est ON.

- 1. Appuyez sur la touche MODE jusqu'à ce que le voyant CON s'allume.
- 2. Appuyez sur la flèche droite jusqu'à ce que "CO2 RLY XXX" apparaisse sur le bloc d'affichage des messages.
- 3. Appuyez sur la flèche vers le haut/bas pour activer (ON) ou désactiver (OFF) cette fonction.
- 4. Appuyez sur ENTER pour enregistrer ce paramètre
- 5. Appuyez sur la touche MODE pour revenir en mode RUN ou sur la flèche vers la droite/gauche pour passer au paramètre suivant/précédent.

Réglage du zéro pour les capteurs CT de CO₂

Lors de l'installation d'un nouveau capteur CT de CO₂, les deux nombres figurant sur l'étiquette collée en usine sur la cellule du capteur doivent être saisis pour étalonner le CO₂ sur l'unité.

Remarque Pour faciliter le travail des techniciens, une étiquette contenant les deux nombres apposés sur la cellule du capteur est fixée à l'intérieur du rack du circuit électronique. ▲

- 1. Appuyez sur la touche MODE jusqu'à ce que le voyant CON s'allume.
- 2. Appuyez sur la flèche droite jusqu'à ce que "T/CZ# XXXX" apparaisse sur le bloc d'affichage des messages.
- 3. Appuyez sur la flèche haut/bas pour remplacer le zéro par le nombre figurant sur l'étiquette.
- 4. Appuyez sur ENTER pour enregistrer ce paramètre.
- 5. Appuyez sur la touche MODE pour revenir en mode RUN ou sur la flèche vers la droite/gauche pour passer au paramètre suivant/précédent.

Réglage du nouvel intervalle de mesure pour les capteurs CT de CO₂

Lors de l'installation d'un nouveau capteur CT de CO₂, les deux nombres figurant sur l'étiquette collée en usine sur la cellule du capteur doivent être saisis pour étalonner le CO₂ sur l'unité.

Remarque Pour faciliter le travail des techniciens, une étiquette contenant les deux nombres apposés sur la cellule du capteur est fixée à l'intérieur du rack du circuit électronique. ▲

- 1. Appuyez sur la touche MODE jusqu'à ce que le voyant CON s'allume.
- 2. Appuyez sur la flèche droite jusqu'à ce que "T/CS# XXXX" apparaisse sur le bloc d'affichage des messages.
- 3. Appuyez sur la flèche haut/bas pour remplacer le nombre affiché par celui figurant sur l'étiquette.
- 4. Appuyez sur ENTER pour enregistrer ce paramètre.
- 5. Appuyez sur la touche MODE pour revenir en mode RUN ou sur la flèche vers la droite/gauche pour passer au paramètre suivant/précédent.

Réglage d'une limite d'alarme basse de HR

Sur les unités où l'option HR est installée, une limite d'alarme basse de HR peut être enregistrée. Cette limite est le pourcentage d'humidité relative à l'intérieur de l'incubateur qui déclenche une alarme basse de HR. Sa valeur peut être définie entre 0 et 90 % d'humidité relative. Le réglage usine est 0 % HR, ce qui désactive l'alarme.

- 1. Appuyez sur la touche MODE jusqu'à ce que le voyant CON s'allume.
- 2. Appuyez sur la flèche droite jusqu'à ce que "RH LO XX" apparaisse sur le bloc d'affichage des messages.
- 3. Appuyez sur la flèche haut/bas pour modifier l'alarme basse de HR.
- 4. Appuyez sur ENTER pour enregistrer la limite d'alarme basse de HR.
- 5. Appuyez sur la touche MODE pour revenir en mode RUN ou sur la flèche vers la droite/gauche pour passer au paramètre suivant/précédent.

Activation des contacts d'alarme de HR

L'alarme basse de HR peut être programmée pour déclencher les contacts d'alarme à distance. La fonction est activée si le paramètre est réglé sur ON. Un réglage sur OFF empêche l'alarme basse de HR de déclencher les contacts. Le réglage usine est ON.

- 1. Appuyez sur la touche MODE jusqu'à ce que le voyant CON s'allume.
- 2. Appuyez sur la flèche droite jusqu'à ce que "RH RLY XXX" apparaisse sur le bloc d'affichage des messages.
- 3. Appuyez sur la flèche vers le haut/bas pour activer (ON) ou désactiver (OFF) cette fonction.
- 4. Appuyez sur ENTER pour enregistrer ce paramètre.
- 5. Appuyez sur la touche MODE pour revenir en mode RUN ou sur la flèche vers la droite/gauche pour passer au paramètre suivant/précédent.

Activation de l'affichage Temp/HR

Sur les unités équipées de l'option HR, le bloc d'affichage supérieur à sept segments du panneau de commande peut être configuré de manière à afficher la température en permanence, à afficher l'humidité relative en permanence ou à basculer de l'une à l'autre. Si l'unité n'est pas équipée de l'option HR, le bloc d'affichage supérieur affiche toujours la température. Si la température est réglée sur ON et l'humidité relative sur OFF, la température est affichée en permanence. Si la température est réglée sur OFF et l'humidité relative sur ON, cette dernière est affichée en permanence. Si les deux paramètres sont réglés sur ON, l'affichage passe de l'une à l'autre. Le réglage usine est par défaut sur ce mode si l'option HR est présente.

- 1. Appuyez sur la touche MODE jusqu'à ce que le voyant CON s'allume.
- 2. Appuyez sur la flèche droite jusqu'à ce que "DISP TMP XXX" ou "DISP RH XXX" apparaisse sur le bloc d'affichage des messages.
- 3. Appuyez sur la flèche vers le haut/bas pour activer (ON) ou désactiver (OFF) cette fonction.
- 4. Appuyez sur ENTER pour enregistrer ce paramètre.
- 5. Appuyez sur la touche MODE pour revenir en mode RUN ou sur la flèche vers la droite/gauche pour passer au paramètre suivant/précédent.



Tableau 3-1. Mode Configuration Page 1/3



Tableau 3-1. Mode Configuration Page 2/3



Tableau 3-1. Mode Configuration Page 3/3

Section 4 Alarmes

Le système d'alarmes des Incubateurs Series 8000DH est répertorié dans le tableau ci-dessous. Lorsqu'une alarme est active, le message s'affiche sur le bloc d'affichage des messages. Une pression sur la touche 🗇 désactive l'alarme sonore pendant l'intervalle de rappel. Néanmoins, l'alarme visuelle continue jusqu'à ce que l'incubateur revienne à une situation normale. Les alarmes sont uniquement momentanées. Lorsqu'une condition d'alarme se produit, puis disparaît, l'incubateur efface automatiquement la condition d'alarme et le message correspondant.

Désignation	Message	Retard	Rappel	Relais
Aucune condition d'alarme	SYSTEM OK ou CLASS 100			
Temp > point de consigne surtempérature	SYS IN OTEMP	0 minutes	15 minutes	Oui
Défaut capteur temp. air (voir Section 4)	AIR SNSR ERR	0 minutes	15 minutes	Non
Panne du contrôleur de température (voir Sect. 4)	TMP CTRL ERR	0 minutes	15 minutes	OUI
Défaut capteur CO2 (voir Section 4)	CO2 SNSR ERR	0 minutes	15 minutes	Non
Période de rappel de changement de filtre HEPA écoulée (voir Section 3)	REPLACE HEPA	0 minutes		Non
Porte intérieure ouverte	DOOR OPEN	15 minutes	15 minutes	Non
CO2 > alarme haute de CO2	CO2 IS HIGH	15 minutes	15 minutes	Programmable
CO2 < alarme basse de CO2	CO2 IS LOW	15 minutes	15 minutes	Programmable
TEMP < alarme basse de TEMP	TEMP IS LOW	15 minutes	15 minutes	Programmable
HR < limite d'alarme basse de HR (option HR)	RH IS LOW	30 minutes	15 minutes	Programmable

Iabicau 4-1. Alainie et description	Tal	bleau	4-1.	Alarme	et	description
-------------------------------------	-----	-------	------	--------	----	-------------

- Tous les retards et intervalles de rappel ont une marge de ±30 secondes -

Lorsque plusieurs conditions d'alarmes se produisent simultanément, les messages actifs s'affichent un par un sur le bloc d'affichage et sont mis à jour toutes les cinq secondes. Une pression sur la touche 🖂 lorsque plusieurs alarmes sont présentes désactive la tonalité de toutes les alarmes, jusqu'au prochain rappel 15 minutes plus tard.

L'alarme basse de température "TEMP IS LOW" est désactivée lorsque le point de consigne de la température est réglé sur 10 °C. Les alarmes de CO₂ sont désactivées lorsque le point de consigne du CO₂ est réglé sur 0,0 %.

Panne du contrôleur de température TMP CNTR ERR

En plus des autres fonctions de sécurité des incubateurs Series 8000DH, un thermostat est inclus pour surveiller la température de l'incubateur. Dans le cas improbable d'une panne du contrôleur de la température, le thermostat éteindrait toutes les résistances à une température de l'incubateur de 160 °C ±5 %. Cette fonction de sécurité est destinée à protéger l'incubateur et non les cultures de cellules ou l'équipement à l'intérieur de la cuve en cas de panne du contrôleur de la température. Si une telle panne se produisait, contactez le service technique ou votre distributeur local.

Alarmes de défauts des capteurs

Le microprocesseur des incubateurs Series 8000DH lit en continu les valeurs de tous les capteurs disponibles afin de garantir qu'ils fonctionnent correctement. S'il détecte une erreur, l'incubateur déclenche une alarme sonore et affiche le message approprié. Dans ce cas, contactez votre distributeur local.

MAINTENANCE PREVENTIVE

Incubateurs

Votre équipement a été minutieusement testé et calibré avant sa livraison. Un entretien préventif régulier est important pour maintenir le fonctionnement correct de votre appareil. L'opérateur doit procéder régulièrement au nettoyage et à l'entretien de routine. Pour des performances et une efficacité maximales, il est recommandé de faire vérifier et calibrer périodiquement l'unité par un technicien de maintenance qualifié. Voici un récapitulatif des tâches d'entretien préventif nécessaires. Reportez-vous à la section concernée du manuel d'utilisation pour plus de détails.

traçables NIST. Pour plus d'informations sur la maintenance préventive ou sur les garanties étendues, veuillez contacter distributor. Thermo Fisher Scientific dispose dans de nombreuses régions de techniciens de maintenance qualifiés utilisant des instruments

La fréquence des nettoyages et des étalonnages dépend de l'utilisation, des conditions environnementales et de la précision requise.

Conseils pour tous les incubateurs :

- N'UTILISEZ PAS d'eau de Javel ou de désinfectants à forte teneur en chlore.
- Utilisez de l'eau <u>stérile</u>, distillée ou déminéralisée.

- Évitez de vaporiser du produit de nettoyage sur le capteur de CO₂.
 - N'utilisez pas de gants talqués pour les cultures de tissus.

États-Unis et Canada 888-213-1790 • Fax : 740-373-4189 • <u>http://www.services.controlenv@thermo.com</u> Millcreek Road, Box 649 • Marietta, Ohio 45750 États-Unis • 740-373-4763

Preventive Maintenance for Series 8000DH Incubators

Refer to Manual Section	Action	Daily	Weekly	Monthly	3 to 6 Months	Yearly
-	Check CO ₂ tank levels.	~				
-	Inspect door latch, hinges and door gasket seal.					~
1	Check water level in the humidity pan, $1/2$ " from top.		>			
2	Verify and document CO ₂ , humidity and temperature calibration, as applicable (See Calibration).					>
	Disinfect the interior of the incubator (See Routine Maintenance).				>	
Ι	Replace HEPA, access port filter, air sample filter, and CO ₂ filters*, if applicable (or as needed).					>
5	Procédez au cycle de décontamination selon les besoins.					
1	Maintenance et inspection de la tuyauterie : inspectez l'ensemble de la tuyauterie et des raccordements et remplacez tout tuyau présentant des signes de dégradation ou de dommages.				>	

Section 5 Entretien courant

Mise en garde Avant d'utiliser une quelconque méthode de nettoyage autre que celles recommandées par le fabricant, l'utilisateur doit s'assurer auprès de ce dernier que la méthode proposée n'endommagera pas l'équipement. ▲

Avertissement Il est de la responsabilité de l'utilisateur de nettoyer immédiatement après tous les déversements accidentels de matières dangereuses. Assurez-vous de suivre les politiques EHS locales en ce qui concerne les équipements de protection individuelle, le nettoyage et l'élimination. ▲

Utilisez un désinfectant approprié, tel que quaternary ammonium. Toutes les pièces et les surfaces doivent être minutieusement nettoyés, rincés à l'eau stérile et séchés.

Avertissement L'alcool, même en solution à 70%, est volatile et inflammable. Ne l'utilisez que dans une zone bien ventilée à l'écart de toute flamme nue. Si un composant est nettoyé à l'alcool, ne l'exposez pas à une flamme nue ou à d'autres risques potentiels. Ne vaporisez pas de solutions inflammables sur le capteur CT. La température interne du capteur de CO₂ atteint environ 150 °C lorsque l'unité est en fonctionnement. Laissez-le refroidir suffisamment avant de le nettoyer. N'utilisez pas de bases fortes ni d'agents caustiques. Bien que résistant à la corrosion, l'inox n'est pas pour autant inaltérable à l'épreuve de la corrosion. N'utilisez pas de solutions d'hypochlorite de sodium (eau de Javel), qui pourraient provoquer des piqûres et de la rouille. ▲

Remarque Si la condensation est visible sur le plafond de la chambre, mais le centre de messagerie n'affiche pas ajouter d'eau, puis le commutateur de niveau de liquide pourrait être défectueux. Appelez les services techniques. ▲

Warning Si l'appareil est en service, débranchez le cordon d'alimentation de l'appareil et la source d'alimentation, éteignez tous les régulateurs de gaz, et laisser l'appareil refroidir avant de désinfecter.

Nettoyage de l'intérieur de l'incubateur

Prévoyez un kit de remplacement du filtre HEPA (voir liste des Pièces de rechange) à portée de main avant de nettoyer l'intérieur de l'incubateur.

Nettoyage de l'intérieur de l'incubateur (cont.)

1. Retirez les étagères, le filtre du port d'accès, le filtre HEPA, le filtre d'échantillonnage de l'air et sa tubulure, ainsi que les panneaux latéraux gauche et droit. Jetez les filtres HEPA et le filtre du port d'accès. Voir Figure 5-1.



Figure 5-1. Emplacements des composants

- 2. Lavez la tubulure du filtre d'échantillonnage de l'air, les étagères et les panneaux latéraux avec un désinfectant, puis rincez-les à l'eau stérile. Vous pouvez également passer les étagères et les panneaux latéraux à l'autoclave.
- 3. Lavez le joint de la porte intérieure avec du désinfectant. Ce joint peut être retiré pour être nettoyé ou remplacé.
- 4. Lavez l'intérieur de l'incubateur avec du désinfectant, en commençant par le haut et en descendant. Reportez-vous aux instructions relatives au désinfectant sur la durée nécessaire avant rinçage. Lavez l'intérieur et l'extérieur de la porte intérieure. La chambre et la porte doivent être rincés à l'eau stérile jusqu'à ce que le désinfectant ait été éliminé. Après le rinçage de la chambre, vaporisez-le à l'alcool à 70 %.
- 5. Installez les panneaux latéraux gauche et droit, le joint de la porte intérieure, le filtre du port d'accès ainsi que celui d'échantillonnage de l'air et sa tubulure et vaporisez-les à l'alcool à 70 %.
- 6. Installez un nouveau filtre HEPA.
- 7. Installez les étagères et vaporisez-les à l'alcool à 70 %.

Nettoyage de l'extérieur de l'incubateur

Nettoyez l'extérieur de l'incubateur à l'aide d'une éponge humide ou d'un chiffon doux bien essoré et d'un détergent doux dissous dans l'eau. Essuyez avec un chiffon doux.

Nettoyage des portes en verre

La porte en verre de la chambre et les portes intérieures indépendantes en verre en option peuvent être nettoyées avec le même désinfectant que celui utilisé à l'intérieur de l'incubateur. Il est impératif de les rincer à l'eau distillée stérile pour éliminer tout résidu de désinfectant. Les portes doivent ensuite être séchées à l'aide d'un chiffon doux.

Certaines précautions doivent être prises lors du nettoyage et de l'entretien des portes en verre de l'incubateur : l'humidité dépose les matériaux alcalins (sodium, Na) à la surface du verre. L'évaporation de l'humidité concentre ces résidus alcalins et peut provoquer des taches ou une opalescence à la surface du verre. Les agents chimiques de nettoyage dont le pH dépasse 9 et la chaleur (autoclave) accélèrent le processus de corrosion. Pour cette raison, il est très important de rincer et d'essuyer les portes de verre après leur nettoyage. Le passage en autoclave des portes de verre doit être évité.

Il n'existe aucune méthode simple pour réparer le verre corrodé. La plupart du temps, le verre doit être remplacé.

Nettoyez le bac humidificateur avec du savon, de l'eau et un désinfectant de laboratoire à usage général. Rincez à l'eau stérile et vaporisez de l'alcool à 70%. Le bac peut être passé à l'autoclave.

Remplacez le filtre HEPA lorsque le rappel REPLACE HEPA s'affiche. Le rappel REPLACE HEPA peut être défini pour déclencher une alarme après une durée spécifiée comprise entre 1 et 12 mois. La durée par défaut du rappel est la durée recommandée de 6 mois. Pour plus de détails, voir Section 3.

Pour accéder au seul fusible de l'incubateur :

- 1. Éteignez l'incubateur avec l'interrupteur et débranchez le cordon secteur.
- 2. Retirez les deux vis en haut du panneau de commande. Saisissez le panneau de commande de chaque côté et tirez-le en ligne droite vers l'extérieur.
- 3. La Figure 5-3 montre l'emplacement du fusible. Voir Tableau 5-1 ciaprès pour connaître les caractéristiques du fusible de remplacement.
- 4. Fixez le panneau de commande et remettez l'unité sous tension. Si le fusible saute lorsque vous aurez remis l'incubateur sous tension, contactez votre distributeur local.

Tableau 5-1. Remplacement des fusibles

Tension du fusible	Réf. fabricant	Courant nominal	Vitesse de rupture	Code lettre CEI
220/240 V	BUSS GMC-0.5A	1250 mA	Retard	Т
220/240 V	BUSS GMC-1.0A	1,0 A	Retard	Т

Nettoyage du bac humidificateur

Maintenance du filtre HEPA

Remplacement des fusibles

Remplacement du filtre d'échantillonnage de l'air

- 1. Branchez une extrémité du filtre d'échantillonnage de l'air à la cannelure du tuyau sur le haut de la cuve.
- 2. Branchez l'autre extrémité du filtre sur la tubulure métallique de l'adaptateur de filtre HEPA (Figure 5-2).



Figure 5-2. Branchement sur la tubulure métallique

Electronique

- 1. Éteignez l'incubateur avec l'interrupteur et débranchez le cordon secteur.
 - 2. Retirez les deux vis à l'arrière du panneau supérieur du boîtier.
 - 3. Faites glisser le panneau supérieur vers l'arrière d'environ deux centimètres et demi et soulevez-le.

Principaux composants

Reportez-vous Figure 5-3. Certains composants représentés sont des options installées à l'usine.



Figure 5-3. Schéma du circuit électronique

Cycle de décontamination

Informations à connaître avant de démarrer un cycle

- Le cycle de décontamination nécessite environ 12 heures : chauffage (2-4 heures), décontamination (2 heures) et refroidissement (6-8 heures). Un temps supplémentaire est nécessaire pour vérifier l'étalonnage de la température et du CO₂ après la fin du cycle.
- Pendant le cycle de décontamination, l'incubateur met à jour la température sur la carte de sortie analogique et sur l'incubateur et fixe le CO₂ à son point de consigne et l'humidité relative à 1 % au-dessus de sa limite basse.
- Un kit de remplacement du filtre HEPA (voir liste des Pièces détachées) doit être placé à portée de main avant le lancement du cycle de décontamination.

Informations sur le cycle

- Un nettoyage préliminaire peut être nécessaire. Pour éviter toutes odeurs et tâches à l'intérieur, sur le matériau étuvé, etc., essuyez toutes les traces d'éclaboussures.
- Une odeur peut apparaître pendant le cycle de décontamination. Il s'agit d'un phénomène normal.
- Le cycle de décontamination n'est pas destiné à stériliser d'autres articles, instruments, etc., du laboratoire.
- Pendant le cycle, la chambre de l'unité devient suffisamment chaude pour faire fondre les échantillons, instruments, récipients, etc. laissés à l'intérieur. Les éléments ci-dessous doivent également être retirés.
 - 1) Filtre HEPA
 - 2) Filtre d'échantillonnage de l'air
 - 3) Filtre du port d'accès
 - 4) Eau du bac humidificateur
 - 5) Sonde d'enregistrement de temp./HR, le cas échéant
 - 6) Capteur IR, le cas échéant

- Filtre Filtre HEPA du port d'accès et sa tubulure avec bouchon Bac humidificateur
- Pendant le cycle de décontamination, une décoloration des matériaux peut se produire. Par exemple, l'inox prend une couleur paille après une certaine période d'exposition à des températures élevées. Ce phénomène est normal.

Points de contrôle

Que faire si ?	Alors
Le cycle ne démarre pas ou se te	ermine prématurément. Vérifiez les alarmes : SYS IN OTEMP, AIR SNSR ERR, TMP CTRL ERR.
Les unités sont empilées	Un cycle de décontamination sur l'une des unités affecte le fonctionnement de l'autre. N'utilisez pas la seconde unité pendant un cycle de décontamination. Le cycle de décontamination peut être effectué simultanément sur les deux unités avec une connexion électrique appropriée.
Aucune action n'est effectuée pe	endant 1 minute après la demande du bloc d'affichage. L'unité revient en mode normal de fonctionnement, SYSTEM OK.
Vous devez interrompre le cycle	en cours. Maintenez le bouton vert de lancement du cycle enfoncé pendant 3 secondes.
L'eau n'est pas retirée du bac hu	midificateur pendant le cycle. De la vapeur est produite et peut provoquer des brûlures.
Les composants répertoriés ne s	ont pas retirés avant le lancement du cycle. Ces composants, qui ne peuvent pas supporter les températures du cycle de décontamination, seront détruits.
L'unité n'est pas mise hors tensi	on avant le retrait du capteur IR, le cas échéant. Le capteur peut être endommagé.
Un capteur IR fictif n'est pas ins	tallé. Le câble du capteur, qui ne peut pas supporter les températures du cycle de décontamination, est détruit.
La porte extérieure est ouverte p	endant les phases de chauffage ou de décontamination. Une alarme de porte extérieure se déclenche : CLOSE DOOR sur le bloc d'affichage, accompagné d'une alarme sonore (qui ne peut être désactivée) et visuelle.
La porte extérieure est ouverte p	endant plus de 20 secondes pendant les phases ci-dessus. Le cycle est annulé et l'unité passe en CANCELED COOL PHASE (phase de refroidissement annulée)*.
La porte extérieure est ouverte pe	endant la phase de refroidissement lorsque la température est supérieure ou égale à 60 °C Une alarme de porte extérieure ouverte se déclenche.
Une coupure de courant se produ	uit pendant la phase de chauffage (HEAT PHASE). La phase HEAT PHASE reprend si la température de la cuve était inférieure à 90 °C lorsque le courant a été coupé. La phase HEAT PHASE reprend si la température était supérieure à 90 °C lors de la coupure de courant et si cette température a baissé de moins de 1 °C. L'unité entre dans une phase d'annulation du refroidissement (CANCELED COOL PHASE)* si la baisse de température dépasse 1 °C.
 * Pour plus d'informations sur la ph 	uit pendant la phase de décontamination (STERILIZATION PHASE). Une phase de chauffage (HEAT PHASE) commence si la température de la cuve n'est pas tombée en dessous de 139 °C. Une phase de décontamination (STERILIZATION PHASE) recommence si la température de la cuve atteint 140 °C. La phase CANCELED COOL PHASE* démarre si la température de la cuve est tombée en dessous de 139 °C. ase CANCELED COOL PHASE voir page 5-12

Thermo Scientific

Cycle de décontamination

Avertissement Le cycle de décontamination chauffe les surfaces intérieures de l'incubateur à 140 °C. Tout contact avec une surface à l'intérieur de la porte externe pendant ce cycle peut provoquer des brûlures. ▲

- 1. Videz l'eau du bac humidificateur et replacez ce dernier à l'intérieur de l'incubateur. Retirez tous les échantillons, instruments, récipients, etc. de la cuve.
- 2. Appuyez sur le gros bouton blanc du cycle de décontamination sur le côté droit de l'unité (Figure 5-4) et maintenez-le enfoncé pendant approximativement 3 secondes jusqu'à ce que le voyant s'allume.



Figure 5-4. Emplacement du bouton de décontamination

- 3. SAISISSEZ le code d'accès le cas échéant. Un code d'accès est recommandé pour éviter un déclenchement accidentel du cycle.
- 4. **Prédécontamination avec capteur de CO**² **TC** Le bloc d'affichage bascule entre "REMOVE HEPAs", "REMOVE WATER" et "PRESS ENTER". Si vous n'appuyez pas sur ENTER dans un délai de 1 minute, le message revient à "SYSTEM OK".
- 5. Retirez le filtre HEPA, le filtre d'échantillonnage de l'air et le filtre du port d'accès (Figure 5-5). La tubulure peut rester à l'intérieur de l'unité pendant le cycle de décontamination. Retirez également la sonde de l'enregistreur temp./HR, le cas échéant. Jetez les filtres.



Figure 5-5. Emplacements des filtres

Prédécontamination avec capteur de CO₂ **IR** - Si un capteur IR est branché, l'affichage bascule entre "POWER OFF" et "REMOVE IR". Si l'unité n'est pas mise hors tension dans un délai de 1 minute, le message revient à "SYSTEM OK". **Remarque** Pour retirer le capteur IR (le cas échéant), commencez par mettre l'unité hors tension. Tournez ensuite les fixations 1/4 de tour sur le conduit supérieur jusqu'à ce que le couvercle soit débloqué. Un capteur fictif est fixé à l'intérieur de la plaque (Figure 5-6). ▲



Figure 5-7. Capteur IR et son câble

6. Saisissez le capteur IR et tirez-le avec précaution vers le bas. Dévissez le câble du capteur IR (Figure 5-6). Mettez le capteur de côté. Pour de plus amples information sur la procédure de désinfection du capteur IR, reportez-vous à l'étape 9. Branchez le capteur fictif sur le câble. Installez le capteur fictif à l'emplacement du capteur IR.

Lorsque l'unité est remise sous tension sans le capteur IR, les messages affichés alternent entre "REMOVE HEPAs", "REMOVE WATER" et "PRESS ENTER". Si vous n'appuyez pas sur "ENTER" dans un délai de 1 minute, les messages alternent entre "POWER OFF" et "REPLACE IR". Lorsque l'unité est remise sous tension avec le capteur IR réinstallé, le message affiché revient à "SYSTEM OK".



Figure 5-8. Messages lorsque le capteur IR est retiré

7. Lorsque la touche ENTER est enfoncée, la phase de chauffage est lancée. Le voyant blanc clignote et le message alterne entre "STERILIZING" et "HEAT PHASE". Pendant cette période, le voyant HEAT est allumé et l'unité se chauffe jusqu'à la température de décontamination.



Figure 5-9. Phase de chauffage

8. Lorsque la température de l'air à l'intérieur de l'unité atteint 140,0 °C, la phase de décontamination commence et le message "STERILIZING" s'affiche.



Figure 5-10. Phase de décontamination

9. Au bout de 2 heures environ, un signal sonore de 5 secondes retentit pour indiquer que la décontamination est terminée. La phase de refroidissement commence. Les messages "STERILIZING" et "COOL PHASE" s'affichent alternativement.



Figure 5-11. Phase de refroidissement

10. Fin du cycle de décontamination avec un capteur CT de CO₂ -Lorsque l'incubateur revient à la température de fonctionnement définie ou à 30 °C si cette dernière est inférieure, les messages alternent entre "CYC COMPLETE", "REPL HEPAs" et "PRESS ENTER". Le voyant blanc reste allumé mais ne clignote plus.

Fin du cycle de décontamination avec un capteur de CO₂ IR -Lorsque l'incubateur revient à la température de fonctionnement définie ou à 30 °C si cette dernière est inférieure, les messages alternent entre "CYC COMPLETE", "POWER OFF" et "REPLACE IR". Le voyant blanc reste allumé mais ne clignote plus.

11. Nettoyez le capteur IR à l'isopropanol ou au désinfectant de surface sans rinçage Lysol. Si vous utilisez l'isopropanol, vaporisez le produit sur le capteur (sans saturation) et laissez sécher. Avec du Lysol, vaporisez sur le capteur (sans saturation) et laissez agir quelques minutes. Essuyez avec chiffon doux propre.

Ne saturez pas le capteur et ne l'immergez pas dans le produit de nettoyage.

Éteignez l'unité, ouvrez la chambre et retirez la plaque du capteur IR. Tirez sur le capteur fictif, dévissez le câble en haut de ce dernier, que vous introduirez ensuite dans la pince de la plaque. Reportez-vous Figure 5-12. Vissez le câble sur le sommet du capteur. Le câble est muni d'un détrompeur et ne peut être inséré que dans un sens. Réinstallez le capteur en le poussant à fond dans l'orifice.

Lorsque l'unité est remise sous tension avec le capteur IR réinstallé, les messages affichés alternent entre "CYC COMPLETE", "REPL HEPAs" et "PRESS ENTER".



Figure 5-12. Capteur IR réinstallé

- 12. Ouvrez la chambre et installez un nouveau filtre HEPA, le filtre d'échantillonnage de l'air et le filtre du port d'accès. Appuyez sur ENTER.
- 13. Lorsque vous appuyez sur ENTER, le voyant s'éteint, le message revient à "SYSTEM OK" et la temporisation de rappel de remplacement du filtre HEPA est réinitialisée.
- 14. Remplissez le bac humidificateur de 3 litres d'eau distillée stérile avec le niveau de pureté recommandé (voir Section 1).
- 15. Laissez la température requise et le niveau de CO₂ se stabiliser au moins 12 heures. Il est recommandé de vérifier régulièrement l'étalonnage de la température et du CO₂ pendant la première semaine de retour en fonctionnement de l'unité.

Mise en garde Les hautes températures du cycle de décontamination peuvent modifier de façon significative la sortie du capteur CT de CO₂. (Ce phénomène est normal et ne reflète pas une détérioration du capteur.) Pour cette raison, il est essentiel de vérifier l'étalonnage du CO₂ avant de remettre l'unité en fonctionnement. ▲

Si vous ne disposez d'aucun instrument indépendant pour vérifier l'étalonnage, vous pouvez suivre la procédure ci-après. Après le cycle de décontamination et avant que le CO₂ ne puisse pénétrer dans la cuve:

- 1) Remplissez et installez le bac humidificateur.
- 2) Laissez l'unité se stabiliser 12 heures à la température de fonctionnement.
- 3) Étalonnez l'écran à 0,0 %.

Remarque Si la phase de refroidissement est annulée, les messages affichés alternent entre "CYC CANCELED" et "COOL PHASE".

- Avec un capteur CT de CO₂ Lorsque l'incubateur revient à la température de fonctionnement définie ou à 30 °C si cette dernière est inférieure, les messages alternent entre "CYC CANCELED", "REPL HEPAs" et "PRESS ENTER". Le voyant blanc reste allumé mais ne clignote plus.
- Avec un capteur de CO₂ IR Lorsque l'incubateur revient à la température de fonctionnement définie ou à 30 °C si cette dernière est inférieure, les messages alternent entre "CYC CANCELED", "POWER OFF" et "REPLACE IR". Le voyant blanc reste allumé mais ne clignote plus. Lorsque l'unité est remise sous tension avec le capteur IR réinstallé, les messages affichés alternent entre "CYC CANCELED", "REPL HEPAs" et "PRESS ENTER".

Section 6 Options installées en usine

Ce chapitre décrit les les options de connexions à des équipement externes installées en usine.

Report d'alarmes

Un ensemble de contacts à relais est fourni pour surveiller les alarmes via une prise de téléphone RJ11 à l'arrière du boîtier. L'emplacement de ce connecteur d'alarmes est représenté Figure 6-5. Un cordon de téléphone 3,6 m (réf. 190388) et un adaptateur RJ11-bornier à vis (réf. 190392) peuvent être commandés auprès de votre distributeur local.

L'alarme distante fournit une sortie NO (normalement ouvert), une sortie NC (normalement fermé) et un contact COM (commun). Reportez-vous Figure 6-1.

Les contacts se déclencher dans des conditions de coupure de courant ou de température excessive. Les contacts peuvent également être programmés pour se déclenchent ou non sur des alarmes de température, de CO₂ et de HR. Reportez-vous Section 3, Mode Configuration.



Figure 6-1. Contacts pour alarme à distance

Affichage de l'humidité

Les incubateurs Series 8000DH sont équipés d'un capteur d'humidité destiné à surveiller l'humidité relative à l'intérieur de la cuve. Ce capteur, qui se trouve dans le panneau supérieur de la cuve, fournit un signal affiché par paliers de 1 % sur le panneau de commande. La lecture de l'humidité peut être affichée en continu ou en alternance avec celle de la température. De plus, une limite d'alarme basse peut être définie pour l'humidité lorsque le bac humidificateur est sec. Reportez-vous à la Section 3, Configuration.

Le capteur peut mesurer l'humidité relative entre 10 % et 100 % avec une précision de ±5 % au dessus de 90 %. Reportez-vous à la Section 2, Étalonnage, pour plus de détails sur l'étalonnage de la mesure d'humidité.

Les niveaux d'humidité dans la cuve peuvent être affectés par les facteurs suivants : niveau d'eau dans le bac humidificateur, fréquence des ouvertures de la porte et position du bac humidificateur au fond de la cuve ou sur une étagère.

Mise en garde Les incubateurs équipés d'un capteur de CO_2 à conductivité thermique ne peuvent mesurer et contrôler avec précision la concentration de CO_2 dans l'incubateur que si le niveau d'humidité relative est constant.

Section 7 Caractéristiques

*Ces caractéristiques techniques sont basées sur une tension nominale de 115 V ou 230 V avec une température ambiante comprise entre 22 °C et 25 °C.

Température

Contrôle
Point de consigneNumérique - Touches à effleurement, 0,1 °C
Gamme+5 °C au-dessus de la température ambiante - 50 °C
Uniformité±0,3 °C @ +37 °C
Alarme de suiviIndicateur (limite basse) programmable par l'utilisateu
Surtempérature
programmable par l'utilisateu
Affichage
Protection Contre la Surchauffe
Type Sécurité, action et indicateur contre les températures extrême
CapteurThermostat, indépendant du système de contrôle de la
températur
Voyant
Humidité Relative
CommandeBac humidificateur - vaporisation naturell
Humidité avec bac
Affichage
AlarmeLimite basse de HR avec moniteur de HR en option
CO_2
Contrôle
CapteurCT ou II
Précision affichage
Gamme
Pression d'arrivée15 PSIG (1 bar
Affichage
Étagères
Dimensions
Construction Acier inoxydable (poli à la bande abrasive sur les deux
côtés
Surface
Maximum par cuve
Loading
Standard
Maximum

Structure

Volume intérieur	
Intérieur	Inox type 304 finition brillante
Extérieur	Acier laminé à froid calibre 18
Joint d'étanchéité de porte ex	térieureSantoprene magnétique, moulé quatre faces
Joint d'étanchéité de porte int	érieure À bourrelet silicone
Isolation	
Connecteurs	
Passage de câble	Bouchon silicone amovible 32 mm
Arrivée CO ₂	Cannelée 6,4 mm
Orifice de prélèvement d'écha	ntillon de gazCannelé, monté à
	l'avant
Caractéristiques électriques	
Modèles 115 volts	
	115 VCA, 50/60 Hz, 1 PH, 9,6 FLA
	(Plage de fonctionnement 90-125 VCA)
Modèles 230 volts	
	230 VCA, 50/60 Hz, 1 PH, 4,4 FLA
	(Plage de fonctionnement 180-250 VCA)
Interrupteur d'alimentation/C	oupe-circuit 2 pôles, 12,0 A
Sortie accessoires	Tension égale à l'entrée du boîtier.
75	watts maximum, courant de fuite 0,5 mA
Contacts pour alarme	Fluctuation de température,
	CO ₂ , coupure de courant, NO et NC
Dimensions de l'appareil	
Extérieures	$\dots .66,3 \text{ cm I x } 100,3 \text{ cm H x } 63,5 \text{ cm P}$
Intérieures	$\dots .54,0 \text{ cm I x } 68,1 \text{ cm H x } 50,8 \text{ cm I'}$
Poids	
Net	
Certifications	
Déclaration de Conformité D	isponible sur Demande

Caractéristiques de sécurité

1
Utilisation à l'intérieur uniquement
AltitudeJusqu'à 2 000 mètres
Température
Humidité80 % HR pour des températures allant jusqu'à 31 °C,
décroissance linéaire jusqu'à 50 % HR à 40 °C sans condensation
Fluctuations de l'alimentation secteur
Les variations de l'alimentation secteur ne doivent pas dépasser ±10 %
de la tension nominale.
Catégorie d'installation II ¹
Degré de pollution 2^2
Classe d'équipement ¹

1 La catégorie d'installation (catégorie de surtension) définit le niveau des surtensions transitoires que l'instrument est conçu pour supporter sans danger. Elle dépend de la nature de la source d'alimentation et de ses moyens de protection contre les surtensions. Par exemple, dans la CAT II qui est la catégorie utilisée pour les instruments dans les installations alimentées par une source d'alimentation comparable à celle des hôpitaux, des laboratoires de recherche et de la plupart des laboratoires industriels, la surtension transitoire attendue est de 2500 V pour une alimentation 230 V et de 1500 V pour une alimentation 120 V.

2 Le degré de pollution définit la quantité de pollution conductrice présente dans le milieu de fonctionnement. Un degré 2 de pollution suppose normalement la seule présence d'une pollution non conductrice telle que de la poussière, à l'exception d'une conductivité occasionnelle produite par la condensation.

Thermo Fisher Scientific 401 Millcreek Road Marietta, Ohio 45750 États-Unis

www.thermofisher.com