

thermoscientific

Vanquish

Compartiments à colonnes VC-C10, VH-C10

Manuel d'utilisation

4827.3201-FR Révision 4.0 • Août 2022





Copyright © 2022 Thermo Fisher Scientific Inc. Tous droits réservés.

Traduction du manuel d'utilisation original

Les descriptions de matériel contenues dans cette révision de manuel correspondent aux appareils : VC-C10-A-03, VH-C10-A-03.

Marques déposées

Acrobat, Adobe et Adobe Reader sont des marques de commerce d'Adobe Systems Incorporated. Microsoft et Windows sont des marques de commerce de Microsoft Corporation. MP35N est une marque de commerce de SPS Technologies. Torx est une marque de commerce de Acument Intellectual Properties, LLC. Toutes les autres marques de commerce citées sont la propriété de Thermo Fisher Scientific Inc. et de ses filiales.

Exclusion de responsabilité

Le présent document accompagne les produits de Thermo Fisher Scientific Inc. lors de l'achat et doit être consulté lors de l'utilisation du produit. Le présent document est protégé par le droit d'auteur ; toute reproduction partielle ou complète de ce document est interdite sans le consentement écrit préalable de Thermo Fisher Scientific Inc.

Le présent manuel a été rédigé en toute conscience. Son contenu peut être modifié à tout moment, sans notification, dans des versions ultérieures.

Thermo Fisher Scientific Inc. ne prétend aucunement que le présent document soit exhaustif, correct et exempt d'erreurs. Thermo Fisher Scientific Inc. n'assume aucune responsabilité pour les erreurs, les défauts, les dommages ou les pertes découlant de l'utilisation du présent document, même si les informations du présent document ont été suivies scrupuleusement.

Le présent document ne fait pas partie du contrat de vente conclu entre Thermo Fisher Scientific Inc. et un client. Le présent document ne régit ni ne modifie aucune condition générale. Si les deux documents présentent des informations contradictoires, les conditions générales prévalent.

Manuel papier uniquement

Imprimé en Allemagne sur du papier ultrablanc 100 % sans chlore, fabriqué selon un procédé écologique ne produisant aucune émission de CO₂.

Adresse du fabricant

Dionex Softron GmbH, Part of Thermo Fisher Scientific, Dornierstrasse 4, D-82110 Germering

Nous contacter

Vous pouvez nous contacter selon plusieurs modalités :

Renseignements concernant les commandes

Pour obtenir des renseignements sur les commandes ou bénéficier du service après-vente pour les produits HPLC, veuillez prendre contact avec le service commercial Thermo Fisher Scientific local. Pour en obtenir les coordonnées, consultez la rubrique Contact Us sur http://www.thermofisher.com.

Assistance technique

Pour obtenir une assistance technique pour les produits HPLC, veuillez prendre contact avec l'assistance technique Thermo Fisher Scientific locale. Pour en obtenir les coordonnées, consultez la rubrique Contact Us sur http://www.thermofisher.com.

Table des matières

1	Uti	lisatio	n du présent manuel	. 9
	1.1	Préser	ntation du présent manuel	10
	1.2	Conve	entions	11
		1.2.1	Conventions des messages de sécurité	11
		1.2.2	Avis spécifiques	11
		1.2.3	Conventions typographiques	12
	1.3	Docur	nentation de référence	13
2	Séc	urité.		15
	2.1	Symbo	oles de sécurité et messages d'avertissement	16
		2.1.1	Symboles de sécurité et messages d'avertissement de ce manuel	16
		2.1.2	Respect des consignes du présent manuel	16
		2.1.3	Symboles de sécurité sur l'instrument	17
		2.1.4	Plaque signalétique	17
	2.2	Utilisa	ition prévue	18
	2.3	Précau	utions de sécurité	19
		2.3.1	Informations générales relatives à la sécurité	19
		2.3.2	Qualification du personnel	20
		2.3.3	Équipements de protection individuelle	20
		2.3.4	Consignes de sécurité électrique	21
		2.3.5	Risques résiduels généraux	22
		2.3.6	En cas d'urgence	24
	2.4	Inform	nations sur les solvants et les additifs	25
		2.4.1	Compatibilité générale	25
		2.4.2	Plages pH autorisées	26
		2.4.3	Concentrations autorisées	26
		2.4.4	Informations complémentaires	26
	2.5	Confo	rmité aux normes et directives	28
3	Pré	senta	tion de l'appareil	29
	3.1	Caract	téristiques du compartiment à colonnes	30
	3.2	Princi	pe de fonctionnement	31
	3.3	Comp	osants internes	33
	3.4	Détec	tion des fuites	34

	3.5	Identif	fication des colonnes (ID de colonne)	35
	3.6	Utilisa	tion	36
4	Déb	allage	2	37
	4.1	Déball	age de l'appareil	38
	4.2	Matér	iel fourni	40
5	Inst	allatio	on	41
	5.1	Consig	nes de sécurité pour l'installation	42
	5.2	Install	ation de l'appareil	43
	5.3	Exigen	ces relatives au lieu d'installation	45
		5.3.1	Alimentation électrique	45
		5.3.2	Cordon d'alimentation	46
		5.3.3	Condensation	46
	5.4	Accès	aux composants internes	47
	5.5	Mise e	en place du matériel	51
		5.5.1	Disposition du système	51
		5.5.2	Fixation du compartiment à colonnes au système	53
		5.5.3	Évacuation des liquides	55
		5.5.4	Connexion de l'appareil	57
		5.5.5	Branchement du cordon d'alimentation	59
		5.5.6	Installation de l'appareil de refroidissement post-colonne (VH-C10 uniquement)	59
		5.5.7	Installation de l'appareil de préchauffage passif	61
		5.5.8	Installation de la vanne de commutation pour colonne	63
		5.5.9	Installation de l'identification des colonnes (ID de colonne)	66
	5.6	Mise e	en place des raccordements fluidiques	68
		5.6.1	Informations et consignes générales	68
		5.6.2	Raccordement des capillaires, des raccords et des tubes	69
		5.6.3	Raccordement et installation de la colonne	71
		5.6.4	Raccordement et installation de l'appareil de préchauffage actif	73
		5.6.5	Raccordement des composants à une vanne de commutation pour colonne	76
	5.7	Mise e	en marche de l'appareil	79
	5.8	3 Configuration de l'appareil dans le logiciel 80		

6	Utilisation				
	6.1	Introduction du chapitre 82			
	6.2	Consignes de sécurité pendant l'utilisation			
	6.3	Éléments de commande			
		6.3.1 Panneau de commande			
		6.3.2	Voyants d'état		
	6.4	Mise s	ous et hors tension		
	6.5	Prépar	ration de l'appareil à l'utilisation		
	6.6	Princip	paux paramètres d'utilisation		
	6.7	Delta d	de température de préparation et durée d'équilibrage		
	6.8	6.8 Optimisation des performances de l'appareil			
		6.8.1	Consignes générales		
		6.8.2	Préchauffage actif		
	6.9	Arrêt d	de l'appareil		
		6.9.1	Arrêt de courte durée (interruption de fonctionnement)		
		6.9.2	Arrêt de longue durée		
		6.9.3	Redémarrage après un arrêt de longue durée		
7	Mai	intena	ance et entretien 101		
	7.4				
	7.1	102 Presentation de la maintenance et de l'entretien			
	7.2	Consignes de sécurité pour la maintenance et l'entretien 10			
	7.3 Règles générales de maintenance et d'entretien				
	7.4 Maintenance de routine et maintenance préventive		enance de routine et maintenance préventive 108		
		7.4.1	Plan de maintenance 108		
		7.4.2	Nettoyage ou décontamination de l'appareil 109		
		7.4.3	Fonctions Predictive Performance 111		
	7.5	Vanne	de commutation pour colonne 112		
		7.5.1	Inspection de la vanne de commutation pour colonne à la recherche de fuites		
		752	Remplacement de la vanne de commutation pour colonne 113		
		7.5.3	Remplacement du capot de la vanne de commutation pour colonne 116		
	7.6	Rempl	acement de l'appareil de refroidissement post-colonne (VH-C10		
		unique	ement) 118		
	7.7	Rempl	acement de l'appareil de préchauffage 120		
		7.7.1	Remplacement de l'appareil de préchauffage actif 120		
		7.7.2	Remplacement de l'appareil de préchauffage passif 122		
	7.8	8 Remplacement des fusibles de l'alimentation principale 123			

	7.9	Remplacement de la porte de l'appareil 125
	7.10	Mise à jour du microprogramme de l'appareil 127
	7.11	Déménagement ou expédition de l'appareil 129
		7.11.1 Préparation de l'appareil au déménagement 129
		7.11.2 Déménagement de l'appareil 130
		7.11.3 Expédition de l'appareil 131
8	Dép	pannage 133
	8.1	Généralités sur le dépannage 134
	8.2	Messages 136
	8.3	Résolution de fuites de liquide 141
9	Car	actéristiques techniques 143
	9.1	Caractéristiques de performance 144
	9.2	Caractéristiques physiques 146
10	Acc	essoires, consommables et pièces de rechange 147
	10.1	Généralités 148
	10.2	Kit d'expédition 149
	10.3	Accessoires en option 150
	10.4	Consommables et pièces de rechange 151
11	Anr	nexe
	11.1	Conformité aux normes et directives 154
		11.1.1 Déclarations de conformité 154
		11.1.2 Conformité à la directive DEEE 155
		11.1.3 Conformité au règlement de la FCC 155
		11.1.4 Historique des versions du manuel 156
	Ind	ex

1 Utilisation du présent manuel

Ce chapitre fournit des informations sur le présent manuel, sur les conventions qui y sont adoptées, ainsi que sur la documentation de référence disponible en complément de ce manuel.

1.1 Présentation du présent manuel

Le présent manuel décrit les caractéristiques techniques et le principe de fonctionnement de votre appareil Vanquish[™]. Il fournit les instructions d'installation, de mise en place, de démarrage, d'arrêt, de fonctionnement, de maintenance et de dépannage.

Il contient également des messages de sécurité, des mentions de mise en garde et des avis spécifiques. Conformez-vous à tous ces messages pour éviter les accidents corporels, les dommages à l'appareil ou la perte de données.

Veuillez noter les points suivants :

- La configuration de l'appareil peut varier ; aussi, toutes les descriptions ne s'appliquent pas nécessairement à votre appareil particulier.
- Si un détail ne s'applique qu'à un modèle (ou à une variante), alors ce dernier est identifié par sa désignation.
- Les illustrations de ce manuel ne sont fournies qu'à des fins de compréhension élémentaire. Elles peuvent varier du modèle de l'appareil ou du composant. Toutefois, cela ne change rien aux descriptions. Aucune réclamation ne peut se fonder sur les illustrations de ce manuel.
- Le compartiment à colonnes est appelé *appareil* ou *compartiment* à *colonnes* dans ce manuel.

1.2 Conventions

Cette section traite des conventions utilisées dans ce manuel.

1.2.1 Conventions des messages de sécurité

Les messages de sécurité et mentions de mise en garde de ce manuel apparaissent comme suit :

- Les messages de sécurité ou mentions de mise en garde qui s'appliquent à l'ensemble du manuel et à toutes ses procédures sont regroupés dans le chapitre Sécurité.
- Les messages de sécurité ou mentions de mise en garde qui s'appliquent à l'ensemble d'une section ou à plusieurs procédures d'une section figurent au début de la section concernée.
- Les messages de sécurité qui ne s'appliquent qu'à une seule section ou procédure figurent dans la section ou procédure concernée. Leur mise en page diffère de celle du texte principal.

Les messages de sécurité sont souvent précédés d'un symbole et/ou d'un mot d'alerte. Le mot d'alerte est composé en majuscules et en gras.

Veillez à bien comprendre et à respecter tous les messages de sécurité figurant dans le présent manuel.

1.2.2 Avis spécifiques

La mise en page des avis spécifiques et des notes indicatives du manuel diffère de celle du texte principal. Ces avis et notes figurent dans des encadrés et sont identifiés au moyen d'un titre. Ce titre est composé en majuscules et en gras.

AVIS

Signale des renseignements jugés nécessaires pour éviter toute détérioration de l'appareil ou tout résultat de test erroné.

NOTE Signale des renseignements d'intérêt général ou des informations utiles pouvant simplifier une tâche ou optimiser les performances de l'appareil.

1.2.3 Conventions typographiques

Les conventions typographiques suivantes s'appliquent aux descriptions de ce manuel :

Entrée et sortie de données

Les éléments suivants apparaissent en gras :

- les entrées effectuées par saisie au moyen du clavier ou par sélection au moyen de la souris ;
- les boutons affichés à l'écran sur lesquels vous cliquez ;
- les commandes saisies au moyen du clavier ;
- les noms, par exemple des boîtes de dialogue, des propriétés et des paramètres.

Dans un souci de concision, les expressions et les chemins d'accès longs sont indiqués selon un format condensé, par exemple : Cliquez sur **Fichier > Enregistrer sous**.

Références et messages

- Les références à la documentation complémentaire apparaissent en *italique*.
- Les messages qui s'affichent à l'écran sont signalés par des guillemets.

Point de vue

Sauf mention contraire, les mots *gauche* et *droite* sont employés dans ce manuel selon le point de vue d'un observateur qui dirigerait son regard vers la façade avant de l'appareil.

Mots très importants

Les mots très importants du texte principal apparaissent en *italique*.

Version électronique du manuel (PDF)

La version électronique (PDF) du manuel comporte de nombreux liens sur lesquels vous pouvez cliquer afin de consulter d'autres sections du manuel. Ces liens incluent :

- les entrées de la table des matières ;
- les entrées de l'index ;
- les renvois (texte en bleu).

1.3 Documentation de référence

D'autres documents de référence sont disponibles en plus du présent manuel d'utilisation.

Documentation sur le matériel

La documentation complémentaire relative au matériel comporte les éléments suivants :

- Manuels d'utilisation des autres modules du système Vanquish
- Manuel d'utilisation du système Vanquish
- Instrument Installation Qualification Operating Instructions

Thermo Fisher Scientific fournit des manuels d'utilisation actualisés sous forme de fichiers PDF (Portable Document Format) auxquels vous pouvez accéder à partir de notre site Web sur les manuels client. Pour ouvrir et lire les fichiers PDF, vous devez disposer d'Adobe™ Reader™ ou Adobe™ Acrobat™.

Naviguez vers le site Web suivant : www.thermofisher.com/ HPLCmanuals

Documentation sur le logiciel

La documentation complémentaire relative au logiciel comporte les éléments suivants :

Aide et documents sur Chromeleon™
 L'Aide de Chromeleon, qui fournit des renseignements exhaustifs, constitue un support de référence complet, quels que soient les aspects du logiciel abordés.

De plus, la documentation suivante est disponible (la disponibilité dépend de la version du logiciel) :

- Guide d'installation
 Pour obtenir des renseignements élémentaires sur l'installation et la configuration de l'appareil, consultez le Guide d'installation.
- Aide de Instrument Configuration Manager
 Pour obtenir des renseignements spécifiques sur un appareil en
 particulier, reportez-vous à l'aide Instrument Configuration
 Manager. Dans Chromeleon 7, les appareils sont appelés
 « modules ».
- Guide de démarrage rapide
 Pour obtenir des renseignements sur les principaux éléments de l'interface utilisateur et une aide étape par étape des procédures les plus importantes, consultez le Guide de démarrage rapide.

• *Carte de référence* Pour obtenir un aperçu concis des procédures les plus importantes, consultez la *Carte de référence*.

NOTE L'*Aide* et les documents sur *Chromeleon* sont fournis avec le logiciel.

Documentation tierce

Vous pouvez également consulter la documentation utilisateur fournie par les autres fabricants de substances et de composants, par exemple les fiches de données de sécurité (FDS).

2 Sécurité

Ce chapitre fournit des consignes de sécurité générales et spécifiques et renseigne sur l'utilisation prévue de l'appareil.

2.1 Symboles de sécurité et messages d'avertissement

2.1.1 Symboles de sécurité et messages d'avertissement de ce manuel

Ce manuel comporte des consignes de sécurité afin d'éviter tout risque de blessure pour les personnes utilisant l'appareil.

Les symboles de sécurité et messages d'avertissement du présent manuel incluent :



Soyez toujours attentif aux consignes de sécurité. N'utilisez pas l'instrument sans avoir compris l'intégralité des consignes de sécurité et réfléchi aux conséquences de vos actions.



ATTENTION

Signale une situation de danger potentiel qui, si elle n'est pas évitée, pourrait entraîner des blessures mineures à modérées.



AVERTISSEMENT

Signale une situation de danger potentiel qui, si elle n'est pas évitée, pourrait entraîner des blessures graves.

2.1.2 Respect des consignes du présent manuel

Respectez les consignes suivantes :

- Lisez ce manuel attentivement avant toute installation ou utilisation de l'appareil, afin de vous familiariser avec l'appareil et ce manuel. Ce manuel comporte des informations importantes relatives à la sécurité des utilisateurs, ainsi qu'à l'utilisation et à la maintenance de l'appareil.
- Gardez toujours le manuel à proximité de l'appareil afin de pouvoir vous y reporter rapidement.
- Conservez ce manuel et transmettez-le à tous les utilisateurs ultérieurs.

Δ

Lisez, comprenez et respectez tous les messages de sécurité et mentions de mise en garde figurant dans le présent manuel.

2.1.3 Symboles de sécurité sur l'instrument

Ce tableau répertorie les symboles de sécurité qui apparaissent sur l'appareil ou sur les étiquettes apposées sur celui-ci. Conformez-vous à toutes les consignes de sécurité présentes dans ce manuel, afin d'éviter tout risque de blessures ou de détérioration de l'appareil.

Symbole	Description
	Indique un danger potentiel. Consultez ce manuel afin d'éviter tout risque d'accident corporel et/ou de dommage à l'appareil.
—	L'appareil est sous tension
0	L'appareil est hors tension
~	Indique un courant alternatif.
	Indique que la surface s'échauffe lors de l'utilisation. Ne touchez pas ces surfaces lorsqu'elles sont chaudes.

2.1.4 Plaque signalétique

La plaque signalétique est placée sur l'appareil à proximité des connexions électriques. Elle indique le numéro de série, le numéro de référence, le nom du module, le numéro de révision (le cas échéant), et la série et le calibre des fusibles.

2.2 Utilisation prévue

L'appareil est destiné à faire partie du système Vanquish.

Le système Vanquish est conçu pour analyser des mélanges de composés dans des solutions d'échantillon.

L'appareil doit être utilisé par une personne qualifiée et dans un environnement de laboratoire uniquement.

L'appareil et le système Vanquish sont réservés exclusivement à la recherche en laboratoire.

Ils ne doivent pas être utilisés à des fins de diagnostic.

Pratiques de laboratoire

Thermo Fisher Scientific recommande au laboratoire dans lequel le système Vanquish est utilisé de suivre les meilleures pratiques pour les analyses LC. Celles-ci comprennent notamment :

- L'utilisation d'étalons appropriés
- L'étalonnage régulier
- La définition et le respect de durées limites de conservation pour tous les produits consommables utilisés avec le système
- L'opération du système conformément au protocole de « test développé en laboratoire », contrôlé et validé par le laboratoire

2.3 Précautions de sécurité

2.3.1 Informations générales relatives à la sécurité

Tous les utilisateurs doivent respecter les consignes générales de sécurité présentées dans cette section, ainsi que tous les messages de sécurité spécifiques et mentions de mise en garde décrits ailleurs dans ce manuel, lors de toutes les phases d'installation, utilisation, dépannage, maintenance, arrêt et transport de l'appareil.



Si l'appareil est utilisé d'une manière non spécifiée par Thermo Fisher Scientific, la protection fournie par l'appareil peut être altérée. Respectez les consignes suivantes :

- Utilisez l'appareil uniquement dans le cadre de ses caractéristiques techniques.
- Employez exclusivement les pièces de rechange, composants supplémentaires, options et périphériques spécifiquement autorisés et certifiés pour l'appareil par Thermo Fisher Scientific.
- Ne réalisez que les procédures décrites dans le présent manuel d'utilisation et dans les documents relatifs à la pompe. Suivez toutes les instructions étape par étape et utilisez les outils recommandés pour la procédure.
- N'ouvrez pas le boîtier de l'appareil et des autres composants, à moins que cela ne soit expressément indiqué dans le présent manuel.
- Thermo Fisher Scientific ne saurait être tenu responsable d'éventuels dommages, matériels ou autres, résultant de l'usage inapproprié ou incorrect de l'appareil. En cas de question concernant l'usage approprié de l'instrument, veuillez contacter Thermo Fisher Scientific avant de poursuivre.

Normes de sécurité

Cet appareil est un instrument appartenant à la classe de sécurité I (équipé d'une borne de mise à la terre). Il a été fabriqué et contrôlé conformément aux normes de sécurité internationales.

2.3.2 Qualification du personnel

Respectez les consignes ci-après, qui portent sur la qualification du personnel chargé de l'installation et/ou de l'utilisation de l'appareil.



Seules des personnes compétentes sont autorisées à mettre en service l'appareil et à effectuer des connexions électriques conformément à la réglementation en vigueur.

- Thermo Fisher Scientific recommande que l'installation soit toujours effectuée par des personnes certifiées par Thermo Fisher Scientific (par souci de concision, désignées par la suite comme « techniciens d'entretien Thermo Fisher Scientific »).
- Si une personne autre qu'un technicien d'entretien agréé par Thermo Fisher Scientific met en service et configure le module, il est de la responsabilité de l'installateur d'assurer la sécurité du module et du système.



Fonctionnement général

L'appareil est destiné à une utilisation dans un environnement de laboratoire par du personnel formé et qualifié.

Tous les utilisateurs doivent connaître les dangers liés à l'appareil et aux substances qu'ils utilisent. Tous les utilisateurs doivent se conformer aux indications figurant sur les fiches de données de sécurité (FDS).

2.3.3 Équipements de protection individuelle

Portez des équipements de protection individuelle et suivez les bonnes pratiques de laboratoire afin de vous protéger contre les substances dangereuses. L'adéquation de l'équipement de protection individuelle dépend des risques présentés par les substances. Pour des conseils concernant les dangers et les équipements de protection requis pour les substances que vous utilisez, reportez-vous à la fiche technique de sécurité et de manipulation des substances fournie par le vendeur.



Une installation permettant de se rincer les yeux ainsi qu'un évier doivent se trouver à proximité de l'appareil. Si une substance, quelle qu'elle soit, entre en contact avec vos yeux ou votre peau, rincez abondamment la zone concernée à l'eau, puis consultez un médecin.

Vêtements de protection

Afin de vous protéger contre les éclaboussures de produits chimiques, les liquides dangereux ou toute autre contamination, portez des vêtements de protection appropriés, telle qu'une blouse de laboratoire.

Équipement de protection oculaire

Afin d'éviter que des projections de liquides n'entrent en contact avec vos yeux, revêtez un équipement de protection des yeux approprié, tel que des lunettes de protection avec écrans latéraux. En cas de risque de projection de liquides, portez des lunettes de sécurité intégrales.

Gants

Afin de vous protéger contre les liquides nocifs et d'éviter de vous blesser lors de la maintenance ou de l'entretien, portez des gants de protection appropriés.

2.3.4 Consignes de sécurité électrique



AVERTISSEMENT—Décharge électrique ou détérioration de l'appareil

Des tensions élevées sont présentes dans l'appareil et peuvent causer des décharges électriques ou endommager l'appareil.

- N'apportez aucune modification aux connexions électriques ou aux bornes de mise à la terre.
- Si vous suspectez la présence d'un dommage électrique quelconque, débranchez le cordon d'alimentation et contactez l'assistance technique de Thermo Fisher Scientific pour obtenir de l'aide.
- N'ouvrez pas le boîtier et ne retirez pas les capots de protection, à moins que cela ne soit expressément indiqué dans le présent manuel.
- Ne placez pas de réservoirs de liquides directement sur l'appareil. Le liquide peut s'écouler dans l'appareil et entrer en contact avec les composants électroniques, provoquant ainsi un court-circuit. Posez plutôt les réservoirs de liquides dans le bac à solvants disponible dans le système Vanquish.

2.3.5 Risques résiduels généraux

Lors de l'utilisation de l'appareil, veuillez prendre en considération les risques résiduels généraux suivants liés au travail avec des substances chimiques :



AVERTISSEMENT—Substances dangereuses

Les solvants, phases mobiles, échantillons et réactifs peuvent contenir des substances toxiques, cancérigènes, mutagènes, infectieuses ou d'autres substances nocives. La manipulation de ces substances peut présenter des risques pour la santé et la sécurité.

- Assurez-vous de bien connaître les propriétés de toutes les substances que vous utilisez. Évitez l'exposition à des substances nocives. Au moindre doute concernant une substance, manipulez-la comme s'il s'agissait d'une substance potentiellement dangereuse.
- Portez les équipements de protection individuelle nécessaires et suivez les bonnes pratiques de laboratoire.
- Réduisez le volume des substances au minimum requis pour l'analyse de l'échantillon.
- Évitez de manipuler les réservoirs de solvant à une hauteur supérieure à la hauteur de la tête.
- N'utilisez pas l'appareil dans un environnement potentiellement inflammable.
- Évitez l'accumulation de substances nocives. Assurez-vous que le lieu d'installation est bien ventilé.
- Éliminez les déchets dangereux de manière écologique, conformément à la réglementation locale en vigueur. Suivez un protocole d'élimination des déchets réglementé et approuvé.



AVERTISSEMENT—Danger biologique

Les matériaux présentant un risque biologique, tels que les microorganismes, les cultures cellulaires, les tissus, les liquides corporels et autres agents biologiques, peuvent transmettre des maladies infectieuses. Pour éviter les infections par ces agents :

- Partez du principe que toutes les substances biologiques sont, tout du moins potentiellement, infectieuses.
- Portez les équipements de protection individuelle nécessaires et suivez les bonnes pratiques de laboratoire.
- Éliminez les déchets biologiques dangereux de manière écologique, conformément à la réglementation locale en vigueur. Suivez un protocole d'élimination des déchets réglementé et approuvé.



AVERTISSEMENT—Auto-inflammation des solvants

Les solvants dont la température d'auto-inflammation est inférieure à 150 °C peuvent s'enflammer au contact d'une surface chaude (par exemple, à la suite d'une fuite dans le système chromatographique).

Évitez d'utiliser ces solvants.



AVERTISSEMENT—Vapeurs dangereuses

Certains échantillons et phases mobiles peuvent contenir des solvants volatils ou inflammables. La manipulation de ces substances peut présenter des risques pour la santé et la sécurité.

- Évitez l'accumulation de ces substances. Assurez-vous que le lieu d'installation est bien ventilé.
- Évitez la proximité des flammes nues et des sources d'étincelles.
- N'utilisez pas l'appareil en présence de gaz ou de fumées inflammables.



ATTENTION—Fuite de substances dangereuses provenant de capillaires en PEEK

Certains capillaires du système sont en PEEK. Le gonflement ou l'exposition à des acides peut provoquer une fuite dans les capillaires en PEEK ou leur éclatement. Certaines substances chimiques, telles que le chloroforme (CHCl₃), le diméthyle sulfoxyde (DMSO) ou le tétrahydrofurane (THF), peuvent provoquer un gonflement du PEEK. Certains acides concentrés, tels que l'acide sulfurique et l'acide nitrique, ou un mélange d'hexane, d'acétate d'éthyle et de méthanol, peuvent endommager le PEEK.

- Ces substances peuvent cependant être utilisées dans le cadre de procédures de nettoyage, à condition que l'exposition soit brève.
- Pour plus d'informations, consultez la documentation technique relative à la résistance chimique du PEEK.



ATTENTION—Réaction allergique

Certains capillaires du système sont composés de MP35N™, un alliage à base de nickel et de cobalt. Le contact avec la peau peut provoquer des réactions allergiques chez les personnes sensibles au nickel ou au cobalt.



ATTENTION—Risque d'étincelles par décharge électrostatique

Les liquides circulant dans les capillaires peuvent générer de l'électricité statique. Cet effet se manifeste en particulier avec les capillaires isolants et les solvants non conducteurs (par exemple, l'acétonitrile pur). Une décharge d'énergie électrostatique peut provoquer la formation d'étincelles, ce qui peut présenter un risque d'incendie.

Évitez de générer de l'électricité statique à proximité du système chromatographique.

2.3.6 En cas d'urgence



AVERTISSEMENT—Risque pour la sécurité

En cas d'urgence, débranchez l'appareil de l'alimentation électrique.

2.4 Informations sur les solvants et les additifs

2.4.1 Compatibilité générale

Afin d'assurer un fonctionnement optimal du système Vanquish, respectez les recommandations suivantes relatives à l'utilisation des solvants et des additifs :

- Le système doit être utilisé exclusivement avec des solvants et additifs compatibles en phase inversée (RP).
- Utilisez uniquement des solvants et additifs compatibles avec tous les éléments du trajet d'écoulement.

NOTE Dans un système Vanquish Core, des solvants et des additifs compatibles avec la phase normale (NP) peuvent être utilisés si les modules du système ont été modifiés pour les applications NP. Consultez le *Manuel d'utilisation du système Vanquish*.

Compatibilité avec le joint de piston

 Dans de rares cas, une durée de vie réduite des joints de piston (UHMW-PE) a été observée en phase inversée après une exposition prolongée à des phases mobiles à pH élevé contenant de l'hydroxyde d'ammonium.

2.4.2 Plages pH autorisées

Système (configuration standard)	Plages pH autorisées	Remarques		
Vanquish Core	1-13	• Valeur de pH de 2 (Vanquish Horizon/Flex) :		
Vanquish Horizon Vanquish Flex	2-12	utilisable à court terme uniquement. La durée d'application doit être aussi brève que possible. Rincez abondamment le système après ces applications.		
		 Valeur de pH de 1-2 (Vanquish Core) : la durée d'application doit être aussi brève que possible. Rincez abondamment le système après ces applications. 		
		 Pour les valeurs de pH supérieures à 9,5 avec des détecteurs optiques : évitez d'utiliser des phases mobiles présentant une valeur de pH supérieure à 9,5 avec des détecteurs optiques. Celles-ci peuvent altérer le fonctionnement et les performances optiques de la cellule à écoulement du détecteur. 		

Plages pH autorisées (configuration standard du système) :

2.4.3 Concentrations autorisées

Concentrations autorisées	(configuration	standard du	système) :
	Comparation	standard du	systemej.

Système (configuration standard)	Chlorure	Tampon	Remarques
Vanquish Core	inférieur ou égal à 0,1 mol/L	inférieur ou égal à 1 mol/L	Avec une concentration élevée en chlorure, la durée d'application doit être aussi brève que
Vanquish Horizon Vanquish Flex	inférieur ou égal à 1 mol/L	-	possible. Rincez abondamment le système après ces applications.

2.4.4 Informations complémentaires

 Pour obtenir des informations détaillées sur les matériaux utilisés dans le trajet d'écoulement analytique de l'appareil, reportez-vous au chapitre *Caractéristiques techniques* de ce manuel. Pour obtenir des informations sur les matériaux utilisés dans le trajet d'écoulement d'autres modules du système Vanquish, consultez le chapitre *Spécifications* dans le *Manuel d'utilisation* du module correspondant.

- Respectez les instructions et recommandations générales sur l'utilisation de solvants et d'additifs dans le système chromatographique. Consultez la section Utilisation de solvants et additifs du Manuel d'utilisation du système Vanquish.
- Consultez également les *Manuels d'utilisation* pour tous les modules du système Vanquish. Ils contiennent des instructions et informations supplémentaires.

AVIS

Si votre configuration de système inclut un détecteur non standard, par exemple un détecteur d'aérosols chargés ou un détecteur d'indice de réfraction, consultez le *Manuel d'utilisation* du détecteur pour des recommandations spécifiques vis-à-vis des solvants et des additifs.

2.5 Conformité aux normes et directives

Thermo Fisher Scientific procède à une évaluation et à des tests complets de ses produits afin de garantir une conformité totale avec les réglementations nationales et internationales en vigueur. Au moment de sa livraison, l'appareil respecte toutes les normes de compatibilité électromagnétique (CEM) et de sécurité applicables, telles que décrites dans le présent manuel.

Toute modification que vous apportez à l'appareil peut potentiellement annuler la conformité avec une ou plusieurs de ces normes de sécurité et de compatibilité électromagnétique. Les modifications de l'appareil comprennent le remplacement d'une pièce ou l'ajout de composants, d'options ou de périphériques non spécifiquement autorisés et certifiés pour le produit par Thermo Fisher Scientific. Afin de garantir une conformité constante avec les normes de sécurité et de CEM, les pièces de rechange et les composants, options et périphériques supplémentaires doivent être commandés auprès de Thermo Fisher Scientific ou de l'un de ses représentants agréés.

L'appareil a été expédié de l'usine dans un état garantissant sa sécurité d'utilisation.

Consultez également

Conformité aux normes et directives () page 154)

3 Présentation de l'appareil

Ce chapitre vous présente l'appareil ainsi que ses principaux composants.

3.1 Caractéristiques du compartiment à colonnes

L'appareil comprend les principales caractéristiques suivantes :

- Thermostatisation des colonnes selon une plage de température comprise entre +5°C et 85°C (VC-C10) ou +5°C et 120°C (VH-C10)
- Refroidissement des colonnes au maximum de 18°C par rapport à la température ambiante, la limite de température inférieure étant +5°C
- Prise en charge des modes de thermostatisation par air calme ou forcé
- Chambre à colonnes pouvant accueillir deux colonnes, avec appareil de préchauffage et précolonne selon les besoins de l'application Longueur maximale possible de l'ensemble : 387 mm
- Identification de colonne, afin d'enregistrer le type de colonne et ses principales propriétés conformément aux bonnes pratiques de laboratoire (BPL)
- Des guides pour capillaires sont situés au-dessus et en-dessous de la chambre à colonnes pour l'acheminement des capillaires
- En option, le compartiment à colonnes peut également comporter :
 - des vannes de commutation pour colonne pour passer rapidement d'une colonne à l'autre ;
 - un appareil de préchauffage pour la thermostatisation de l'éluant avant sa pénétration dans la colonne ;
 - un appareil de préchauffage actif avec une véritable régulation de la température pour chauffer l'éluant avant sa pénétration dans la colonne à une température définie par l'utilisateur
- Appareil de refroidissement post-colonne (VH-C10 uniquement) Pour refroidir l'éluat avant qu'il n'entre dans le détecteur

Pour connaître les caractéristiques techniques, reportez-vous à Caractéristiques de performance (> page 144).

3.2 Principe de fonctionnement

Le critère fondamental d'un compartiment à colonnes destiné à des applications de chromatographie en phase liquide est la capacité à maintenir la température de consigne avec le plus de précision possible. Cependant, pour obtenir des résultats de mesure de bonne qualité, la stabilité de la température est plus importante que la précision de la température. Grâce à ses circuits électroniques de pointe, ce compartiment à colonnes permet de stabiliser la température selon une précision de ±0,1°C.

La conception à optimisation thermique du compartiment à colonnes réduit la durée d'équilibrage de la température entre la colonne et l'éluant.

Les éléments thermoélectriques réchauffent ou refroidissent les composants situés à l'intérieur de la chambre à colonnes. Un appareil de préchauffage actif ou passif peut être installé pour porter l'éluant à la température de la colonne avant que l'éluant ne pénètre dans celle-ci. Cela permet d'éviter les gradients de température dans la première partie de la colonne et une dégradation des performances de séparation.

VH-C10 uniquement :

De plus, un appareil de refroidissement post-colonne peut être installé afin de refroidir l'éluat avant qu'il n'entre dans le détecteur. Le refroidissement post-colonne minimise le bruit du détecteur et optimise les résultats de la détection.

Modes de thermostatisation

Le compartiment à colonnes prend en charge les modes de thermostatisation de colonnes suivants :

- Mode air calme
 - En mode air calme, la colonne se réchauffe sur toute sa longueur.
 - La température à l'intérieur de la colonne est identique à celle de l'extérieur de la colonne.
 - L'éluant et l'échantillon se déplacent à la même vitesse dans la colonne qu'à l'extérieur de celle-ci (absence d'accélération du débit).

Le mode air calme limite les fronts diffus et les traînées de pic.

- Mode air forcé
 - En mode air forcé, la surface de la colonne est maintenue à la température de consigne sur toute la longueur de la colonne.
 - La température à l'intérieur de la colonne est supérieure à celle de l'extérieur de la colonne.
 - L'éluant et l'échantillon se déplacent plus rapidement dans la colonne qu'à l'extérieur de celle-ci (accélération du débit).

Il en résulte des pics plus larges et une dégradation possible des performances de séparation.

Les illustrations représentent la distribution thermique à l'intérieur de la colonne pour les deux modes :



Illustration 1: Conditions thermiques à l'intérieur de la colonne en mode air calme



Illustration 2: Conditions thermiques à l'intérieur de la colonne en mode air forcé

N°	Description
1	Direction du flux dans la colonne
2	Profil du flux
3	Température constante sur la surface de la colonne
4	Augmentation de la température dans la colonne selon la direction du flux

Le déplacement de l'éluant dans la colonne induit un frottement dans la colonne dont l'intensité augmente avec le débit. Ce frottement provoque un échauffement de la colonne. Cet effet est particulièrement important avec les débits élevés des applications UHPLC. Plus le débit est élevé, plus le frottement généré est important et plus la colonne s'échauffe.

3.3 Composants internes

Les composants du compartiment à colonnes accessibles par l'utilisateur se situent directement derrière la porte avant :





N°	Description
1	Ports de connexion pour connecteurs d'ID de colonne
2	Vanne de commutation pour colonne supérieure (si installée)
3	VH-C10 uniquement : appareil de refroidissement post-colonne (si installé)
4	Guides pour capillaires
5	Chambre à colonnes (avec et sans capot de chambre)
6	Emplacements de montage d'appareils de préchauffage passifs
7	Ports de connexion pour appareil de préchauffage actif
8	Vanne de commutation pour colonne inférieure (si installée)
9	Bac de fuite avec capteur pour détecter les fuites de liquide
10	Panneau de commande doté de quatre boutons, pour le pilotage de l'instrument

3.4 Détection des fuites

Les fuites constituent un risque potentiel pour la sécurité.

Le capteur de fuites situé à l'intérieur de l'appareil détecte toute fuite de liquide issue des raccordements fluidiques. Le liquide est recueilli dans le bac de fuite prévu à cet effet, puis canalisé vers l'orifice d'évacuation. Cet orifice permet d'évacuer le liquide par le dispositif d'évacuation du système Vanquish.

Lorsque le capteur de fuites détecte une fuite, les voyants d'état deviennent rouges et un signal sonore se déclenche afin de vous alerter. Suivez les instructions de ce manuel pour trouver et éliminer la source de la fuite.

3.5 Identification des colonnes (ID de colonne)

Le système d'identification des colonnes (ID de colonne) vous permet de stocker des informations spécifiques sur un connecteur d'ID de colonne. Lorsque la colonne est installée dans la chambre et que l'identification de colonne est active, les informations de colonne sont mises à jour en continu et peuvent être consultées par l'utilisateur à partir du système de gestion de données chromatographiques.

Pour une vue d'ensemble de l'état de la colonne conforme aux bonnes pratiques de laboratoire, les informations suivantes sont disponibles :

Propriétés primaires de colonne

Les propriétés primaires de colonne permettent une identification fiable de la colonne. Elles incluent, par exemple, l'identifiant de produit, le numéro de lot, le numéro de série et la date de fabrication.

• Propriétés secondaires de colonne

Les propriétés secondaires de colonne fournissent des informations complémentaires, par exemple le fabricant, le remplissage, la granulométrie, les températures minimale et maximale autorisées ou encore les pressions minimale et maximale autorisées.

• Propriétés opérationnelles de colonne

Les propriétés opérationnelles de colonne fournissent des informations sur le cycle de vie de la colonne, notamment le nombre total d'injections exécutées sur la colonne, la date de la première injection et la date d'injection la plus récente. Le système de gestion de données chromatographiques détermine et met à jour les informations opérationnelles à chaque injection. Les propriétés opérationnelles de colonne sont en lecture seule.

3.6 Utilisation

L'appareil a été conçu pour être piloté par un ordinateur équipé du système de gestion de données chromatographiques Chromeleon. Le logiciel Chromeleon assure de façon complète le pilotage de l'instrument ainsi que l'acquisition et la gestion des données.

Pour obtenir une description élémentaire du pilotage de l'instrument et de l'analyse automatisée des échantillons avec le logiciel Chromeleon, consultez le *Manuel d'utilisation du système Vanquish*. L'*Aide de Chromeleon* comporte des renseignements détaillés sur le pilotage et le fonctionnement de la pompe.

NOTE L'appareil peut également être piloté à partir d'autres systèmes de gestion de données, tels que Thermo Scientific[™] Xcalibur[™]. Dans ce cas, il est nécessaire d'installer d'autres logiciels en plus du logiciel de gestion de données. Pour obtenir des renseignements complémentaires, veuillez prendre contact avec le service commercial de Thermo Fisher Scientific.
4 Déballage

Ce chapitre comporte des renseignements sur le déballage de l'appareil, ainsi que sur les éléments fournis.

4.1 Déballage de l'appareil

Emballage endommagé, constatation d'un défaut à la réception

Inspectez la boîte de transport afin d'y déceler d'éventuels signes de détérioration externe puis, une fois le produit déballé, inspectez l'appareil afin d'y déceler d'éventuels signes d'une détérioration mécanique qui aurait pu survenir pendant le transport.

Si vous soupçonnez que l'appareil a été endommagé pendant le transport, informez-en immédiatement le transporteur et Thermo Fisher Scientific. L'assurance transport n'est valide que si la détérioration est immédiatement signalée.

Déballage

AVIS

Pour lever ou transporter l'appareil, saisissez-le par les côtés. Ne déplacez et ne soulevez jamais l'appareil par la porte avant. Cela risque d'endommager les portes ou l'appareil.

NOTE Conservez la boîte de transport ainsi que l'ensemble du matériel d'emballage. Vous en aurez besoin si vous déménagez ou expédiez l'appareil.

Outils nécessaires

Tournevis type Torx[™] T20

Procédez comme suit

- 1. Placez la boîte de transport par terre, puis ouvrez-la.
- 2. Retirez le kit d'expédition de la boîte.
- 3. Saisissez l'appareil par les côtés.
- 4. Sortez l'appareil de sa boîte de transport en le soulevant doucement et délicatement.
- 5. Placez l'appareil sur une surface stable, sur son panneau latéral gauche, avec le fond de l'appareil dirigé vers vous.
- Le cas échéant
 Retirez tout emballage supplémentaire. Laissez tous les films
 protecteurs recouvrant l'appareil en place jusqu'à ce qu'il soit
 correctement positionné dans le système empilé.

 Repérez les 2 pieds et les 4 vis de fixation dans le kit de transport. Le pied comportant l'orifice d'évacuation est le pied avant de l'appareil.

NOTE Vous pouvez fixer le pied avec l'orifice d'évacuation dirigé vers l'avant ou l'arrière de l'instrument.

8. À l'aide du tournevis, serrez les 2 vis pour monter le pied avant.



Illustration 4: Montage des pieds

- 9. À l'aide du tournevis, serrez les 2 vis pour monter le pied arrière.
- 10. Saisissez les côtés supérieur et inférieur de l'appareil, puis soulevez l'instrument afin de le placer en position verticale.
- 11. Transportez l'appareil jusqu'à l'emplacement où il doit être installé.
- 12. Certaines surfaces, dont la porte de l'appareil, sont recouvertes d'un film protecteur lors du transport. Retirez les films protecteurs de toutes les surfaces recouvertes.



ATTENTION—Renversement

Le compartiment à colonnes peut se renverser.

Pour éviter toute blessure corporelle ou détérioration de l'appareil, fixez systématiquement le compartiment à colonnes au boîtier du système Vanquish. Pour obtenir des instructions détaillées, reportez-vous à la section Fixation du compartiment à colonnes au système (> page 53).

4.2 Matériel fourni

Le matériel suivant est fourni :

- Compartiment à colonnes
- Kit d'expédition
- Manuel d'utilisation (téléchargeable à partir du site Web sur les manuels client)
- Cordon d'alimentation

Pour obtenir des informations concernant le kit d'expédition ou le réapprovisionnement en pièces, reportez-vous à la section Accessoires, consommables et pièces de rechange (▶ page 147).

5 Installation

Ce chapitre précise les exigences relatives au lieu d'installation et décrit comment mettre en place, installer et configurer l'appareil dans le système Vanquish et dans le logiciel de chromatographie.

5.1 Consignes de sécurité pour l'installation

Veuillez respecter les consignes de sécurité suivantes :



Respectez tous les messages d'avertissement et mentions de mise en garde figurant à la section Précautions de sécurité (> page 19).



ATTENTION—Renversement

Le compartiment à colonnes peut se renverser.

Pour éviter toute blessure corporelle ou détérioration de l'appareil, fixez systématiquement le compartiment à colonnes au boîtier du système Vanquish. Pour obtenir des instructions détaillées, reportez-vous à la section Fixation du compartiment à colonnes au système (> page 53).



ATTENTION—Décharge électrique ou détérioration de l'appareil

Lorsque l'appareil est éteint, il n'est pas complètement isolé électriquement tant que le cordon d'alimentation est branché. Réaliser des réparations alors que l'appareil est connecté à l'alimentation électrique peut entraîner des accidents corporels.

- Débranchez toujours le cordon d'alimentation avant de procéder à des travaux de réparation à l'intérieur de l'appareil.
- Si vous devez enlever des capots ou des panneaux, ne branchez pas le cordon d'alimentation à l'appareil tant que les panneaux et les capots sont démontés.

AVIS

Pour lever ou transporter l'appareil, saisissez-le par les côtés. Ne déplacez et ne soulevez jamais l'appareil par la porte avant. Cela risque d'endommager les portes ou l'appareil.

5.2 Installation de l'appareil

Un technicien d'entretien agréé par Thermo Fisher Scientific installe et configure le système Vanquish, notamment tous les modules et toutes les options ou pièces qui les accompagnent. Le technicien d'entretien vérifie que l'installation est correcte et que le système et les modules Vanquish fonctionnent conformément aux spécifications. Le technicien fait également une démonstration du fonctionnement de base et des principales fonctionnalités du système.

Si une personne autre qu'un technicien d'entretien agréé par Thermo Fisher Scientific installe l'appareil, il convient de procéder comme suit.

AVIS

L'appareil fait partie du système Vanquish Aussi, suivez la séquence d'installation des modules du système exposée dans le *Manuel d'utilisation du système Vanquish*.

- Suivez les consignes de sécurité et respectez l'ensemble des exigences relatives au lieu d'installation. Reportez-vous aux sections Consignes de sécurité pour l'installation (▶ page 42) et Exigences relatives au lieu d'installation (▶ page 45).
- Mettez en place le matériel de l'appareil. Reportez-vous à la section Mise en place du matériel (▶ page 51).
- 3. Mettez en place les raccordements fluidiques. Consultez la section Mise en place des raccordements fluidiques (▶ page 68).
- 4. Mettez l'appareil sous tension. Reportez-vous à la section Mise en marche de l'appareil (▶ page 79).

NOTE

Avant de mettre un module du système Vanquish sous tension pour la première fois, vérifiez que le logiciel de chromatographie est installé sur l'ordinateur du système de données. Les pilotes USB nécessaires sont recherchés automatiquement à la mise sous tension et le système d'exploitation Windows™ peut détecter l'appareil.

5. Configurez l'appareil dans le logiciel. Reportez-vous à la section Configuration de l'appareil dans le logiciel (▶ page 80).

 Recommandation : Procédez à une qualification de l'installation de l'instrument.

Sous Chromeleon, un assistant vous guide dans la procédure de qualification. Sur la **Chromeleon 7 Console** : Cliquez sur **Tools > Instrument Qualification > Installation Qualification**.

Suivez les instructions du manuel *Instruments Installation Qualification Operating Instructions*. Ce manuel fournit des informations sur les éléments matériels nécessaires ainsi que des instructions détaillées.

AVIS

Si l'appareil est utilisé avec un autre système de gestion de données, consultez la documentation du logiciel que vous utilisez et/ou exécutez la qualification manuellement. Le manuel *Instruments Installation Qualification Operating Instructions* fournit des informations sur les paramètres à adapter et sur les paramètres nécessaires.

 Recommandé : procédez à une qualification opérationnelle. Le kit de qualification comporte l'ensemble des éléments matériels nécessaires à la qualification ainsi que des instructions détaillées.

Déplacement de l'appareil après l'installation

Si vous devez déplacer l'appareil après sa mise en place et son installation dans le système Vanquish, préparez-le en vue du transport, puis déménagez-la. Suivez les instructions sous Déménagement ou expédition de l'appareil (▶ page 129).

5.3 Exigences relatives au lieu d'installation

Un environnement d'utilisation approprié est indispensable pour assurer une performance optimale de l'appareil.

Cette section présente les principales exigences concernant le lieu d'installation. Veuillez noter les points suivants :

- Utilisez l'appareil uniquement dans des conditions de laboratoire appropriées.
- L'appareil est destiné à faire partie du système Vanquish. Respectez les exigences relatives au lieu d'installation pour le système Vanquish telles qu'indiquées dans le Manuel d'utilisation du système Vanquish.
- Pour obtenir les caractéristiques techniques, reportez-vous à la section Caractéristiques techniques (> page 143) et consultez la section Caractéristiques techniques des Manuels d'utilisation pour les autres modules du système Vanquish.
- Pour connaître les risques résiduels, consultez la section Risques résiduels généraux (> page 22).

5.3.1 Alimentation électrique

L'appareil s'adapte à un vaste éventail d'alimentations électriques ; il accepte toutes les tensions secteur de la plage spécifiée pour l'appareil.



ATTENTION—Décharge électrique ou détérioration de l'appareil

Le branchement de l'appareil à une tension secteur supérieure ou inférieure à celle spécifiée présente un risque de blessure corporelle pour l'utilisateur ou de détérioration pour l'appareil.

Branchez l'appareil uniquement à la tension spécifiée.

5.3.2 Cordon d'alimentation

Les cordons d'alimentation sont conçus pour respecter les spécifications des prises murales du pays dans lequel ils sont utilisés. L'extrémité du cordon d'alimentation branchée dans la prise électrique de l'appareil est identique pour tous les cordons d'alimentation. L'extrémité du cordon d'alimentation qui est branchée dans la prise murale est différente.



AVERTISSEMENT—Décharge électrique ou détérioration de l'appareil

- N'utilisez jamais de cordons d'alimentation autres que ceux fournis par Thermo Fisher Scientific pour l'appareil.
- Utilisez uniquement un cordon d'alimentation conçu pour le pays dans lequel vous utilisez l'appareil.
- N'utilisez pas de rallonges électriques.
- Ne branchez jamais l'appareil à une prise électrique partagée avec d'autres instruments (prises multiples par exemple).
- Utilisez l'appareil uniquement avec une prise d'alimentation avec borne de terre protectrice.
- En cas d'urgence, le cordon d'alimentation doit être facilement accessible à tout instant afin de pouvoir débrancher l'appareil de l'alimentation électrique.



AVERTISSEMENT—Décharge électrique ou détérioration d'un produit

Une mauvaise utilisation des cordons d'alimentation peut entraîner des blessures corporelles ou une détérioration de l'appareil. Utilisez les cordons d'alimentation fournis par Thermo Fisher Scientific uniquement pour l'usage auquel ils sont destinés. Ne les utilisez pas à d'autres fins, par exemple pour connecter d'autres instruments.

5.3.3 Condensation

AVIS—La présence de condensation dans l'appareil peut endommager les composants électroniques.

- Prévenez ou minimisez les conditions pouvant entraîner l'accumulation de condensation dans l'appareil lors de son utilisation, de son expédition ou de son stockage. Par exemple, évitez les changements rapides ou importants des conditions de l'environnement de travail.
- Si vous suspectez la présence de condensation, laissez l'appareil se réchauffer à température ambiante. Cela peut prendre plusieurs heures. Patientez jusqu'à disparition complète de la condensation avant de brancher l'appareil à l'alimentation électrique.

5.4 Accès aux composants internes

Pour accéder aux composants internes du compartiment à colonnes, ouvrez la porte du compartiment à colonnes, puis retirez le capot de la chambre à colonnes.

Ouverture et fermeture de la porte du compartiment à colonnes

- 1. Pour ouvrir la porte, tirez légèrement dessus.
- 2. Pour fermer la porte, poussez dessus.

Ouverture de la chambre à colonnes



ATTENTION—Surfaces chaudes

Les surfaces internes de l'appareil peuvent s'échauffer lors de l'opération. Ne touchez pas les surfaces chaudes pour éviter de vous brûler.

- Ne touchez pas la face interne du capot de la chambre à colonnes. Lorsque vous ouvrez le capot de la chambre à colonnes, saisissez-le uniquement par le porte-capot.
- Attendez que les surfaces chaudes aient refroidi avant de commencer toute procédure de remplacement ou de maintenance.
- 1. Ouvrez la porte du compartiment à colonnes.
- Tournez les deux verrous de 180° dans le sens antihoraire, jusqu'à ce qu'ils se bloquent en position verticale. Un déclic vous confirme que les verrous sont en position correcte pour l'ouverture de la chambre à colonnes.



Illustration 5: Déverrouillage du capot de la chambre à colonnes

- 3. Saisissez le capot de la chambre à colonnes par le porte-capot.
- 4. Tirez légèrement le capot de la chambre à colonnes vers vous puis poussez-le vers la droite avec la face interne du capot orientée vers la face interne de la porte.



Illustration 6: Ouverture du capot de la chambre à colonnes

Fermeture de la chambre à colonnes

- 1. Saisissez le capot de la chambre à colonnes par le porte-capot.
- 2. Tirez légèrement le capot de la chambre à colonnes vers vous puis poussez-le vers la gauche avec la face interne du capot orientée vers la chambre à colonnes.



Illustration 7: Fermeture du capot de la chambre à colonnes

3. Avant de fermer la chambre à colonnes, alignez soigneusement les 4 paumelles du capot de la chambre à colonnes pour permettre sa rotation.



Illustration 8: Paumelles du capot de la chambre à colonnes

4. Tournez les deux verrous de 180° dans le sens horaire, jusqu'à ce qu'ils se bloquent en position verticale.
Repérez la mention Closed située au-dessus des verrous, et veillez à ce que les verrous soient alignés sur ce repère.

NOTE Utilisez toujours le compartiment à colonnes avec la porte du compartiment à colonnes et le capot de chambre à colonnes fermés. Dans le cas contraire, vous ne pourrez pas démarrer des analyses.

5.5 Mise en place du matériel

Cette section, qui traite de la mise en place du matériel, comporte des renseignements sur les raccords, les connecteurs et les câbles de l'appareil.

5.5.1 Disposition du système

L'appareil fait partie du système Vanquish. Les modules sont généralement placés dans le système empilé, leur disposition dépendant de la configuration du système.



Illustration 9: Système Vanquish, configuration standard (exemple)

N°	Description
1	Bac à solvants
2	Détecteur
3	Passeur d'échantillon
4	Pompe
5	Socle du système
6	Compartiment à colonnes

Dans la configuration standard, le système Vanquish comporte un seul compartiment à colonnes, installé sur le côté droit du système.

Pour obtenir des instructions sur la mise en place du système empilé, consultez le *Manuel d'utilisation du système Vanquish*.

NOTE Une configuration étendue ou une configuration avec le compartiment à colonnes installé sur le côté gauche est également disponible. Veuillez noter ce qui suit :

- La configuration étendue sur le côté droit peut inclure jusqu'à trois compartiments à colonnes. Jusqu'à deux kits de conversion sont requis pour assembler la configuration étendue à plusieurs compartiments à colonnes.
- Le compartiment à colonnes peut également être installé sur le côté gauche du système empilé Vanquish. Dans ce cas, l'assemblage du compartiment à colonnes nécessite un kit de conversion. Chaque compartiment à colonnes supplémentaire installé sur le côté gauche nécessite un kit de conversion supplémentaire. Lorsqu'un chargeur Vanquish ou que le Vanquish Display sont utilisés avec le système Vanquish, le compartiment à colonnes doit être installé sur le côté droit du système.
- Pour obtenir des renseignements sur les commandes, veuillez prendre contact avec le service commercial Thermo Fisher Scientific local.



Installation par un technicien d'entretien uniquement

Seuls les techniciens d'entretien certifiés par Thermo Fisher Scientific sont habilités à installer la configuration étendue comprenant jusqu'à trois compartiments à colonnes ou à installer le compartiment à colonnes sur le côté gauche du système.

5.5.2 Fixation du compartiment à colonnes au système

NOTE Avant de fixer au système le compartiment à colonnes, raccordez la conduite d'évacuation à l'orifice d'évacuation du système. Pour obtenir des instructions, consultez le *Manuel d'utilisation du système Vanquish*.

Éléments nécessaires

- 4 écrous rectangulaires en T
- 2 pièces transversales
- 4 vis
- 4 entretoises

Outils nécessaires

Tournevis type Torx T20

Procédez comme suit

1. Alignez les trous de vis des deux écrous rectangulaires en T, des entretoises et de la pièce transversale.



Illustration 10: Barre de montage

N°	Description
1	Entretoise
2	Encoche
3	Vis
4	Écrou rectangulaire en T
5	Pièce transversale

- 2. À l'aide des vis, fixez à la pièce transversale les écrous rectangulaires en T et les entretoises. Les écrous rectangulaires en T doivent être positionnés face à face, comme indiqué ci-dessus.
- 3. Faites glisser les écrous rectangulaires en T de la barre de montage dans le rail du boîtier du système, puis poussez-les jusqu'en butée. Les encoches situées sur les barres de montage doivent être dirigées vers le haut. Assurez-vous que la pièce transversale est orientée comme indiqué dans l'illustration.

0

4. Serrez les vis sur les barres de montage.





N°	Description
1	Rail
2	Barre de montage

- 5. Soulevez le compartiment à colonnes afin de le placer en position verticale.
- 6. Pour insérer les boutons de maintien dans les encoches de la barre de montage, saisissez le compartiment à colonnes par ses côtés, puis soulevez-le légèrement.



Illustration 12: Fixation du compartiment à colonnes au système empilé

5.5.3 Évacuation des liquides

Éléments nécessaires

- Tube d'évacuation
- Coude
- Bac d'évacuation suivant les besoins
- Coupe-tube (en option)

Procédez comme suit

1. Raccordez le tube d'évacuation au coude.

 Raccordez le coude à l'orifice d'évacuation.
 L'orifice d'évacuation est situé dans la partie avant basse de l'appareil, au-dessous de la porte avant.



Illustration 13: Raccordement du tube d'évacuation

N°	Description
1	Orifice d'évacuation
2	Coude

 Placez l'extrémité libre du tube d'évacuation dans un bac d'évacuation adéquat.

Si vous devez raccourcir le tube, utilisez un coupe-tube. Veillez à ce que la coupe soit à angle droit par rapport à l'axe du tube et que l'extrémité ne présente aucune irrégularité.

AVIS

Pour maintenir une évacuation appropriée, placez le bac d'évacuation plus bas que le compartiment à colonnes. Dans le cas contraire, le liquide pourrait refluer à l'intérieur de l'instrument et l'endommager.

Pour obtenir des renseignements sur l'évacuation des liquides dans le dispositif d'évacuation Vanquish, consultez le *Manuel d'utilisation du système Vanquish*.

5.5.4 Connexion de l'appareil

Connecteurs de l'appareil

Les connecteurs suivants sont fournis sur l'appareil:



Illustration 14: Connecteurs

N°	Description
1	Plaque signalétique indiquant le numéro de série, le numéro de référence, le nom du module, le numéro de révision (le cas échéant), la série et le calibre des fusibles
2	Interrupteur d'alimentation principal (marche/arrêt)
3	Porte-fusible
4	Connecteur d'entrée d'alimentation
5	USB Port USB (Universal Serial Bus) (connecteur de type « B ») Permet de connecter la pompe à d'autres modules du système Vanquish ou à l'ordinateur sur lequel est installé le système de gestion des données, tel que le logiciel Chromeleon
6	Port System Interlink Permet de commander la mise sous tension et hors tension de l'appareil depuis le socle du système Vanquish et la communication de l'appareil (par exemple pour le Vanquish User Interface).

NOTE Thermo Fisher Scientific recommande d'utiliser les ports USB uniquement comme décrit ci-dessus. Si les ports USB sont utilisés à toute autre fin, Thermo Fisher Scientific n'est pas en mesure de garantir un fonctionnement correct.

Procédez comme suit

AVIS

- N'utilisez jamais de câbles de communication défectueux. Si vous suspectez qu'un câble est défectueux, remplacez-le.
- Afin d'assurer un fonctionnement sans accroc, utilisez uniquement les câbles fournis par Thermo Fisher Scientific pour connecter l'appareil.
- 1. Placez l'appareil dans le système tel que requis dans la configuration du système. Pour obtenir plus de détails, consultez le *Manuel d'utilisation du système Vanquish*.
- 2. Connectez les câbles d'interface requis à l'appareil. Pour obtenir des renseignements complémentaires sur le branchement de l'appareil à d'autres modules dans le système Vanquish ou à l'ordinateur comportant le système de gestion des données de chromatographie, consultez le *Manuel d'utilisation du système Vanquish*.
- 3. Branchez le cordon d'alimentation (consultez la section Branchement du cordon d'alimentation (▶ page 59)).

5.5.5 Branchement du cordon d'alimentation

AVIS

La présence de condensation dans un appareil peut endommager les composants électroniques.

- Assurez-vous de l'absence de condensation dans les appareils avant de les brancher à l'alimentation électrique.
- Si vous suspectez la présence de condensation, laissez l'appareil se réchauffer à température ambiante. Patientez jusqu'à disparition complète de la condensation avant de poursuivre.
- 1. Vérifiez que l'interrupteur d'alimentation principal de l'appareil est positionné sur OFF (hors tension).
- 2. Branchez le cordon d'alimentation au connecteur d'entrée d'alimentation sur l'appareil.
- 3. Branchez l'extrémité libre du cordon d'alimentation à une source d'alimentation appropriée.

5.5.6 Installation de l'appareil de refroidissement post-colonne (VH-C10 uniquement)

Pour réduire la température du liquide à la sortie la colonne, installez un appareil de refroidissement post-colonne. Le refroidissement postcolonne minimise le bruit du détecteur et optimise les résultats de la détection.

AVIS

Sur les systèmes Vanquish dotés d'un détecteur muni d'une cellule à écoulement LightPipe™, utilisez toujours un appareil de refroidissement post-colonne afin d'abaisser la température du liquide à la sortie de la colonne à un niveau inférieur à 50 °C. Des températures plus élevées détériorent la cellule à écoulement.

Éléments nécessaires

Appareil de refroidissement post-colonne

Procédez comme suit

- 1. Ouvrez la porte du compartiment à colonnes.
- Fixez l'appareil de refroidissement post-colonne à la plaque de montage située au-dessus de la chambre à colonnes.
 Vous pouvez installer l'appareil de refroidissement post-colonne sur l'un ou l'autre côté, selon l'emplacement du détecteur :
 - si le détecteur est situé à gauche du compartiment à colonnes, le capillaire fixé sur le côté de l'appareil de refroidissement postcolonne doit être dirigé vers la gauche ;
 - si le détecteur est situé à droite du compartiment à colonnes, le capillaire de l'appareil de refroidissement post-colonne doit être dirigé vers la droite ;
 - vous pouvez retourner l'appareil de refroidissement postcolonne : si l'éluat chaud sortant de la colonne vient de la vanne de commutation pour colonne supérieure, le capillaire doit être dirigé vers le haut, et si l'éluat chaud vient directement de la colonne, le capillaire doit être dirigé vers le bas. Si l'éluat chaud de la colonne provient de la colonne, le capillaire doit être dirigé vers le bas.



Illustration 15: Montage d'un appareil de refroidissement post-colonne

- 3. Pour fixer l'appareil de refroidissement post-colonne à la plaque de montage, serrez les 2 vis moletées.
- Configurez l'appareil de refroidissement post-colonne dans le système de gestion des données chromatographiques : Dans la boîte de dialogue Instrument Configuration Manager pour le compartiment à colonnes, activez l'appareil de refroidissement postcolonne.

NOTE Lorsque vous enlevez l'appareil de refroidissement post-colonne, désactivez-le dans la boîte de dialogue Instrument Configuration Manager pour le compartiment à colonnes.

5.5.7 Installation de l'appareil de préchauffage passif

Éléments nécessaires

Appareil de préchauffage passif

Outils nécessaires

Tournevis type Torx T10

Procédez comme suit

 Placez l'appareil de préchauffage passif à l'emplacement désiré. Quatre emplacements différents sont disponibles dans la chambre à colonnes.

NOTE L'appareil de préchauffage passif peut être installé dans n'importe lequel des quatre emplacements illustrés ci-dessous. Il peut être orienté selon les besoins, avec le capillaire de sortie dirigé vers le haut ou le bas, selon la position de l'entrée de la colonne.





N°	Description
1	Emplacements pour appareil de préchauffage passif dans la loge de colonne de gauche
2	Emplacements pour appareil de préchauffage passif dans la loge de colonne de droite

2. Alignez les trous de vis sur l'appareil de préchauffage passif avec ceux de l'emplacement de montage.

3. À l'aide du tournevis, serrez les deux vis et rondelles pour monter l'appareil de préchauffage passif à l'intérieur de la chambre à colonnes.





N°	Description
1	Appareil de préchauffage passif (installé)
2	Vis et rondelles de l'appareil de préchauffage passif

- 4. Raccordez le capillaire de sortie de l'appareil de préchauffage passif à l'entrée de la colonne.
- 5. Raccordez le capillaire d'entrée de l'appareil de préchauffage passif au passeur d'échantillon ou à la vanne de commutation pour colonne.

NOTE L'appareil de préchauffage passif n'est pas contrôlé par le système de gestion de données chromatographiques. Il n'est donc pas nécessaire de configurer cet appareil dans le système de gestion de données chromatographiques. L'appareil de préchauffage passif peut être utilisé pour porter l'éluent à la température de la colonne avant que l'éluant ne pénètre dans celle-ci.

5.5.8 Installation de la vanne de commutation pour colonne

Pour un basculement rapide entre les colonnes, le compartiment à colonnes peut accueillir une ou deux vannes de commutation pour colonne. Ces vannes peuvent être installées aux emplacements indiqués dans la section Composants internes (> page 33).

Éléments nécessaires

Vanne de commutation pour colonne

Outils nécessaires

Tournevis type Torx T10

Préparatifs

- 1. Mettez le compartiment à colonnes hors tension. S'il est sous tension, la vanne de commutation peut être endommagée pendant l'installation.
- 2. Ne retirez pas encore les bouchons de raccordement qui couvrent les orifices des vannes.

Retrait du capot de la vanne de commutation pour colonne

- 1. Ouvrez la porte du compartiment à colonnes.
- 2. À l'aide du tournevis, desserrez les 4 vis du capot de la vanne de commutation.



Illustration 18: Vis du capot de la vanne de commutation

3. Retirez les vis et conservez-les pour une installation ultérieure.

- 4. Retirez le capot et le joint d'étanchéité de la vanne de commutation.



N°	Description
1	Joint
2	Capot

5. Détachez avec précaution le câble de l'arrière du joint d'étanchéité pour vanne de commutation.

NOTE Conservez le capot et le joint d'étanchéité de la vanne de commutation pour couvrir l'ouverture de la vanne de commutation si elle est à nouveau retirée.

Installation de la vanne de commutation pour colonne

- 1. Contrôlez l'orientation du câble. La broche d'alignement située au milieu du connecteur du câble doit être dirigée vers le bas.
- 2. Branchez le câble au connecteur situé à l'arrière de la vanne de commutation pour colonne.
- 3. Enfoncez le câble dans le boîtier.

4. Enfoncez délicatement la vanne de commutation pour colonne dans le boîtier.



Illustration 20: Installation de la vanne de commutation pour colonne

- 5. À l'aide du tournevis, serrez les 4 vis de la vanne de commutation pour colonne.
- 6. Retirez les bouchons de raccordement qui couvrent les orifices des vannes.
- Connectez les conduites de liquide, selon les besoins de l'application. Pour les instructions, reportez-vous à la section Raccordement des composants à une vanne de commutation pour colonne (▶ page 76).
- 8. Mettez le compartiment à colonnes sous tension.
- 9. Configurez la vanne de commutation pour colonne dans le système de gestion de données chromatographiques :
 - a) Dans la boîte de dialogue Instrument Configuration Manager, sélectionnez l'emplacement (**Upper Valve**, **Lower Valve**).
 - b) Sélectionnez le modèle de vanne.
- 10. Mettez la vanne dans les différentes positions, puis testez les raccordements fluidiques.

5.5.9 Installation de l'identification des colonnes (ID de colonne)

L'identification de colonne (ID de colonne) permet de créer le cycle de vie d'une colonne (historique de colonne). Le compartiment à colonnes prend en charge l'identification simultanée de quatre colonnes. Les ports de l'ID de colonne sont situés en regard de la vanne de commutation pour colonne supérieure.

Éléments nécessaires

- Connecteur d'ID de colonne
- Rivet





N°	Description
1	Connecteur d'ID de colonne
2	Rivet
3	Ruban

Procédez comme suit

- 1. Ouvrez la porte du compartiment à colonnes.
- 2. Mettez en place le ruban sur la colonne :
 - a) Enroulez le ruban autour de la colonne.
 - b) Faites passer la tige du rivet à travers un orifice, puis faites pression afin de joindre les deux pièces.



Illustration 22: Ruban avec rivet

3. Branchez le connecteur d'ID de colonne au port d'ID de colonne correspondant.



Illustration 23: Branchement du connecteur d'ID de colonne au port d'ID de colonne

Activation de l'identification de colonnes dans Instrument Configuration Manager

- Lors de la première configuration du compartiment à colonnes, sur la page General, pour le compartiment à colonnes, cochez la case Use Column ID system.
- 2. Sur la page **Columns**, activez les connecteurs d'ID de colonne respectifs.

5.6 Mise en place des raccordements fluidiques

Cette section décrit la mise en place des raccordements fluidiques depuis et vers l'appareil, et si nécessaire, de raccordements fluidiques supplémentaires.

5.6.1 Informations et consignes générales

Conformez-vous aux règles et recommandations suivantes lors de la mise en place des raccordements fluidiques :



Les raccordements fluidiques peuvent contenir des substances dangereuses. Respectez les messages d'avertissement et mentions de mise en garde figurant à la section Précautions de sécurité (> page 19).

- Les composants sales peuvent contaminer le système chromatographique. La contamination diminue les performances des modules et du système dans son ensemble, ou peut même causer une détérioration des modules et du système. Par conséquent :
 - Portez toujours des gants appropriés.
 - Placez les composants uniquement sur une surface propre et non pelucheuse.
 - Veillez à maintenir vos outils propres.
 - Utilisez uniquement des chiffons non pelucheux pour le nettoyage.
- Pour obtenir des instructions et consignes d'installation ainsi que des recommandations relatives à la manipulation, reportez-vous à la section Raccordement des capillaires, des raccords et des tubes (> page 69).

AVIS

Lorsque vous installez des appareils ou des composants dans le système, rincez-les et évacuez le liquide de rinçage avant de les raccorder au trajet d'écoulement du système. Pour rincer les modules Vanquish, suivez les instructions du *Manuel d'utilisation du système Vanquish*.

Procédez comme suit

Pour mettre en place les raccordements fluidiques et terminer l'installation du compartiment à colonnes, procédez comme suit :

- Raccordez puis installez les colonnes et l'appareil de préchauffage (reportez-vous à la section Raccordement et installation de l'appareil de préchauffage actif (> page 73)).
- Raccordez les composants à la vanne de commutation pour colonne (reportez-vous à la section Raccordement des composants à une vanne de commutation pour colonne (▶ page 76)).

5.6.2 Raccordement des capillaires, des raccords et des tubes

Cette section traite de la connexion et de la manipulation des capillaires, des raccords et des tubes.

5.6.2.1 Consignes générales

Lorsque vous raccordez les capillaires et les tubes, suivez les recommandations générales suivantes :

- N'utilisez que les capillaires et les tubes (par exemple, les conduites de solvant ou la tubulure d'évacuation) fournis avec le produit ou bien des capillaires et tubes de complément ou de rechange recommandés par Thermo Fisher Scientific.
- Les raccords doivent être débarrassés de tout contaminant. Même des particules infimes peuvent endommager le système ou fausser les résultats des tests.
- N'installez pas des capillaires ou des tubes ayant subi une contrainte, entaillés, déformés ou présentant toute autre détérioration.
- Installez les capillaires et raccords uniquement aux emplacements pour lesquels ils sont prévus.

5.6.2.2 Raccordement des capillaires Viper

Cette section traite de la procédure de raccordement des capillaires Viper[™]. Tous les raccordements fluidiques Viper du système Vanquish ont été conçus pour être serrés à la main.

Pour connecter les capillaires Viper munis d'une molette, procédez comme suit :

AVIS

- Serrez ou desserrez les capillaires Viper *uniquement* à la main.
 N'utilisez aucun outil autre que la molette fournie avec le capillaire.
- Afin d'éviter d'endommager le capillaire ou le raccord, serrez et desserrez les capillaires Viper *uniquement* lorsque la pression du système est redescendue à zéro.



Illustration 24: Raccord Viper avec molette

N°	Description
1	Molette
2	Capillaire
3	Fente

- 1. Insérez le capillaire Viper dans l'orifice de raccordement.
- 2. Serrez le raccord au moyen de la molette.

NOTE Notez bien la fente pratiquée dans la molette. Vous pouvez facilement retirer la molette des capillaires au moyen de cette fente lorsque l'espace est restreint.

3. Contrôlez l'absence de fuite du raccord. Si le raccord fuit, suivez les étapes ci-dessous. Résolution des fuites des raccords Viper avec molettes

- 1. Serrez davantage le raccord.
- 2. Si la fuite persiste, retirez le capillaire.
- 3. Nettoyez soigneusement les extrémités du capillaire au moyen d'une serviette en papier non pelucheuse humectée d'isopropanol.
- 4. Réinstallez le capillaire.
- 5. Si la fuite persiste, installez un nouveau capillaire Viper.

5.6.3 Raccordement et installation de la colonne

À l'intérieur de la chambre à colonnes, une cloison centrale sépare les loges de colonne gauche et droite. Utilisez un porte-colonne pour fixer chaque colonne à la cloison centrale située à l'intérieur de la chambre à colonnes.



Illustration 25: Porte-colonnes

N°	Description
1	Porte-colonne permettant de placer la colonne à gauche
2	Porte-colonne permettant de placer deux colonnes à la même hauteur
3	Porte-colonne permettant de placer la colonne à droite

Procédez comme suit

1. Raccordez les capillaires à l'entrée et à la sortie de la colonne.

AVIS

Pour éviter toute fuite, vérifiez que tous les raccordements d'entrée et de sortie de la colonne sont correctement effectués lors de la pose ou du remplacement des capillaires.



2. Fixez le porte-colonne au raccord de la colonne.

Illustration 26: Fixation du porte-colonne au raccord de la colonne

3. Fixez le porte-colonne sur la cloison centrale située à l'intérieur de la chambre à colonnes.

Le porte-colonne peut être fixé n'importe où sur la cloison centrale.



Illustration 27: Mise en place de la colonne à l'intérieur de la chambre à colonnes

AVIS

Pour éviter d'endommager la cloison centrale, ne faites pas coulisser les supports le long de la cloison. Si vous souhaitez modifier leur position, retirez-les et repositionnez-les comme souhaité.
4. Acheminez les capillaires par les guides pour capillaires. Ces derniers sont situés au-dessus et en dessous de la chambre à colonnes.



Illustration 28: Acheminer le capillaire par le guide pour capillaires

5.6.4 Raccordement et installation de l'appareil de préchauffage actif

Cette section traite du raccordement et de la mise en place de l'appareil de préchauffage actif dans le compartiment à colonnes.

Éléments nécessaires

- Porte-colonne
- Appareil de préchauffage actif
- Filtre en ligne Viper (facultatif)

Préparatifs

AVIS

L'appareil de préchauffage actif est rempli d'isopropanol lorsque le compartiment à colonnes est expédié depuis le site de fabrication. Un isopropanol de qualité insuffisante peut endommager la colonne.

- Rincez les composants remplis d'isopropanol au moyen de solvants miscibles à l'isopropanol. À défaut, utilisez un solvant intermédiaire de polarité similaire.
- Installez un raccord-union Viper dans le trajet d'écoulement pour remplacer la colonne pendant la procédure de rinçage.

Procédez comme suit

NOTE Utilisez un filtre en ligne Viper entre l'appareil de préchauffage actif et la colonne afin d'éviter que des particules non dissoutes ne pénètrent dans la colonne.

- 1. Raccordez le filtre en ligne Viper à la colonne.
- 2. Raccordez l'appareil de préchauffage actif au filtre en ligne Viper.
- 3. Fixez le porte-colonne au raccord de l'appareil de préchauffage actif.
- 4. Fixez le porte-colonne sur la cloison centrale située à l'intérieur de la chambre à colonnes.



Illustration 29: Mise en place de l'appareil de préchauffage et de la colonne à l'intérieur de la chambre à colonnes

- 5. Avant de connecter pour la première fois le câble de préchauffage au port de l'appareil de préchauffage, vérifiez que l'appareil de préchauffage est bien activé dans la boîte de dialogue Instrument Configuration Manager du compartiment à colonnes. Une fois l'appareil de préchauffage activé dans le système de gestion de données chromatographiques, il détecte automatiquement la connexion de l'appareil de préchauffage.
- Connectez le câble de préchauffage au port de l'appareil de préchauffage.
 Les ports de connexion d'appareils de préchauffage sont situés à côté de la vanne de commutation pour colonne inférieure.

NOTE Utilisez le port de connexion gauche si l'appareil de préchauffage est installé dans la loge de colonne gauche et le port droit s'il est installé dans la loge droite.



Illustration 30: Port de connexion de l'appareil de préchauffage

7. Passez le câble de l'appareil de préchauffage par le serre-câble situé juste au-dessus de l'appareil de préchauffage. Chaque câble préchauffage et guide-câble se trouve au-dessus du port de préchauffage. Si le câble de préchauffage est trop long, passez le câble dans le guide.





N°	Description
1	Guide-câble
2	Câble de l'appareil de préchauffage

5.6.5 Raccordement des composants à une vanne de commutation pour colonne

Plusieurs types de vannes de commutation pour colonne sont disponibles pour le compartiment à colonnes. Pour obtenir les informations de commande, reportez-vous à la section Accessoires en option (▶ page 150). Lors de leur expédition, les vannes de commutation pour colonne sont accompagnées d'informations détaillées expliquant les différents raccordements fluidiques possibles.

L'installation d'une vanne de commutation pour colonnes est décrite à la section Installation de la vanne de commutation pour colonne (> page 63).

Cette section présente l'exemple des raccordements fluidiques pour une vanne de commutation pour colonne à 2 positions et 6 ports dans le compartiment à colonnes Vanquish.

Éléments nécessaires

- Vanne de commutation pour colonne
- Capillaires Viper, selon les besoins de l'application
- Appareil de refroidissement post-colonne (facultatif)
- Appareil de préchauffage actif ou passif (facultatif)
- Capuchons Viper

AVIS

Du liquide peut se déverser si des ports de la vanne restent ouverts. Le capteur de fuites du compartiment à colonne détecte tout déversement de liquide et signale une fuite au moyen d'un message d'erreur et d'une alarme sonore.

Pour éviter tout déversement, bouchez systématiquement les ports de vanne non utilisés à l'aide de capuchons Viper.

5.6.5.1 Raccordements fluidiques pour l'utilisation d'une vanne de commutation pour colonne

La figure présente les raccordements fluidiques recommandés pour l'utilisation d'une vanne de commutation pour colonne.



Illustration 32: Exemple : raccordements sur une vanne de commutation pour colonne 2 pos. 6 ports. Emplacement de vanne supérieur. Gauche : position 6-1 ; droite : position 1-2

Procédez comme suit

- 1. Raccordez le capillaire Viper provenant de la pompe ou du passeur d'échantillon au port 6 de la vanne de commutation pour colonne.
- Selon la position de la colonne, raccordez un appareil de préchauffage actif ou passif au port 1 ou 5 de la vanne de commutation pour colonne et raccordez l'appareil de préchauffage à l'entrée de la colonne.
- 3. Raccordez un capillaire Viper au port 2 ou 4 de la vanne de commutation pour colonne et à la sortie de la colonne.
- Raccordez un appareil de refroidissement post-colonne ou un capillaire Viper au port 3 de la vanne de commutation pour colonne et à l'entrée de la cellule à écoulement du détecteur.

5.6.5.2 Raccordements fluidiques pour l'utilisation de deux vannes de commutation pour colonne

La figure présente les raccordements fluidiques recommandés pour l'utilisation de deux vannes de commutation pour colonne.



Illustration 33: Exemple : mise en place de deux vannes de commutation pour colonne dans le trajet d'écoulement. Gauche : position 6-1 ; droite : position 1-2

Procédez comme suit

- 1. Raccordez le capillaire Viper provenant de la pompe ou du passeur d'échantillon au port 3 de la vanne de commutation pour colonne inférieure.
- Selon la position de la colonne, raccordez un appareil de préchauffage actif ou passif au port 2 ou 4 de la vanne de commutation pour colonne inférieure et raccordez l'appareil de préchauffage à l'entrée de la colonne.
- 3. Raccordez un capillaire Viper au port 1 ou 5 de la vanne de commutation pour colonne supérieure et à la sortie de la colonne.
- Raccordez un appareil de refroidissement post-colonne ou un capillaire Viper au port 6 de la vanne de commutation pour colonne supérieure et à l'entrée de la cellule à écoulement du détecteur.

5.7 Mise en marche de l'appareil

NOTE

Avant de mettre un module du système Vanquish sous tension pour la première fois, vérifiez que le logiciel de chromatographie est installé sur l'ordinateur du système de données. Les pilotes USB nécessaires sont recherchés automatiquement à la mise sous tension et le système d'exploitation Windows™ peut détecter l'appareil.

Pour mettre en marche l'appareil, procédez comme suit :

- Vérifiez que le bouton d'alimentation situé sur la partie avant gauche du socle du système Vanquish (bouton d'alimentation du système) est enfoncé. Si le bouton d'alimentation n'est pas enfoncé (proéminent), appuyez dessus pour mettre sous tension le socle du système.
- 2. Mettez l'appareil sous tension au moyen de son interrupteur d'alimentation principal.

Mettez l'appareil hors tension au moyen de l'interrupteur d'alimentation principal lorsque vous en recevez l'instruction, par exemple pendant des opérations de maintenance. Il ne suffit pas d'appuyer sur le bouton d'alimentation du système pour couper complètement l'alimentation électrique de l'appareil.

Consultez également

B Mise sous et hors tension (▶ page 87)

5.8 Configuration de l'appareil dans le logiciel

Dans ce manuel, il est admis que le logiciel de chromatographie est déjà installé sur l'ordinateur du système de données et qu'une licence en cours de validité y est disponible.

Pour obtenir des renseignements complémentaires sur la configuration du système Vanquish dans le logiciel, reportez-vous au *Manuel d'utilisation du système Vanquish*.

L'Aide du logiciel que vous utilisez comporte des renseignements détaillés sur les paramètres de chaque page de propriétés.

6 Utilisation

Ce chapitre décrit les éléments nécessaires au pilotage de l'appareil et vous informe sur les opérations de routine et sur l'arrêt de l'appareil.

6.1 Introduction du chapitre

Dans ce chapitre, il est admis que la configuration initiale de l'appareil a déjà été effectuée. Dans le cas contraire, reportez-vous aux instructions de la section Installation () page 41).

Pour obtenir une description élémentaire du pilotage de l'instrument et de l'analyse automatisée des échantillons avec le logiciel Chromeleon, consultez le *Manuel d'utilisation du système Vanquish*. L'*Aide de Chromeleon* comporte des renseignements détaillés sur le pilotage et le fonctionnement de la pompe.

Les descriptions logicielles de se manuel se rapportent à Chromeleon 7. La terminologie peut différer de celle utilisée dans les autres versions logicielles.

6.2 Consignes de sécurité pendant l'utilisation

Lorsque vous utilisez l'appareil veuillez respecter les consignes de sécurité suivantes :



Respectez tous les messages d'avertissement et mentions de mise en garde figurant à la section Précautions de sécurité (> page 19).



ATTENTION—Surfaces chaudes

Les surfaces internes de l'appareil peuvent s'échauffer lors de l'opération. Ne touchez pas les surfaces chaudes pour éviter de vous brûler.

- Attendez que les surfaces chaudes aient refroidi avant de les toucher.
- Pour refroidir rapidement les surfaces chaudes situées à l'intérieur du compartiment à colonnes, appuyez sur le bouton FAST COOL du panneau de commande. Pour en savoir plus, reportez-vous à la section Panneau de commande (> page 84).

AVIS

Prenez aussi en considération les consignes de sécurité suivantes :

- Pour éviter des dommages causés par des fuites ou d'avoir utilisé la pompe à sec, réglez toujours la limite de pression inférieure de la pompe.
- Si une fuite ou les signes d'une fuite sont détectés, arrêtez le débit de pompe et remédiez immédiatement au problème.

6.3 Éléments de commande

L'appareil a été conçu pour être piloté principalement depuis un ordinateur équipé du logiciel de chromatographie.

De plus, les éléments suivants sont également disponibles sur l'appareil :

- Panneau de commande
 Les boutons du panneau de commande vous permettent d'exécuter certaines fonctions directement depuis l'appareil.
- Voyants d'état
 Les LED (diodes électroluminescentes) de la barre d'état, située sur la façade avant de l'appareil, et la LED STATUS, situées sur le panneau de commande, permettent un contrôle visuel rapide de l'état opérationnel de l'appareil.

6.3.1 Panneau de commande

Le panneau de commande, situé à l'intérieur de l'appareil, vous permet d'exécuter certaines fonctions directement depuis celui-ci. Lorsque vous appuyez sur un bouton, un bref signal sonore confirme que la fonction s'exécute.

Lorsque l'appareil est connecté sous le logiciel Chromeleon, il se peut que certaines fonctions ne soient pas disponibles à partir du panneau de commande (reportez-vous aux renseignements ci-dessous, dans la présente section).



Illustration 34: Panneau de commande

STATUS

La LED **STATUS** permet un contrôle visuel rapide de l'état opérationnel de l'appareil.

Lorsque la porte est fermée, la barre à LED située sur la façade avant indique l'état opérationnel.

Pour obtenir plus de détails, reportez-vous à la section Voyants d'état (> page 86).

MUTE ALARM		
	Un signal sonore vous prévient lorsque l'appareil détecte un problème, par exemple une fuite. Pour désactiver le signal sonore de l'alarme en cours, appuyez sur ce bouton. Éliminez la source de l'alarme en 10 minutes. Sinon, le signal sonore reprend. Si l'appareil détecte un problème différent, le signal sonore vous prévient de nouveau immédiatement.	
FAST COOL		
	Ce bouton permet de refroidir rapidement les surfaces chaudes du compartiment à colonnes, par exemple pour permettre un accès rapide et sans danger aux composants afin de les remplacer ou effectuer une opération de maintenance. Pour obtenir plus de détails, reportez-vous à la section Principaux paramètres d'utilisation () page 90).	
	Pendant l'opération de refroidissement, la LED du bouton est rouge. Une fois la température de consigne atteinte, la LED devient verte, ce qui indique que vous pouvez accéder aux composants en toute sécurité.	
	Pour annuler le mode Fast Cool, appuyez de nouveau sur le bouton. La LED du bouton s'éteint et les composants sont chauffés pour être portés à la dernière température de consigne. Le mode Fast Cool peut également être annulé via le logiciel Chromeleon.	
UPPER VALVE , LOW	/ER VALVE	
	Ces boutons permettent de faire passer les vannes de commutation pour colonne (supérieure ou inférieure) d'une position à l'autre (si installées).	
	Les LED situées à côté des boutons indiquent la position actuelle des vannes à 2 positions et 6 ports. La LED supérieure est verte lorsque la vanne est en position 1 et la LED inférieure est verte lorsque la vanne est en position 2. Les LED sont éteintes pendant la commutation de la vanne.	
Lorsque l'appareil est connecté sous le logiciel Chromeleon		
	Lorsque l'appareil est connecté sous le logiciel Chromeleon, le fonctionnement des boutons est comme suit :	

- Aucune injection ou séquence n'est en cours : Tous les boutons du panneau de commande sont actifs.
- Une injection ou séquence est en cours : Seul le bouton MUTE ALARM reste actif depuis le panneau de commande, ce qui vous permet de désactiver le signal sonore de l'alarme en cours.

6.3.2 Voyants d'état

La barre à LED d'état, située sur la façade avant de l'appareil, et la LED **STATUS**, située sur le panneau de commande intérieur, fournissent des informations sur l'état de l'appareil.

Barre à LED

Les couleurs de la barre à LED donnent les informations suivantes :

Barre à LED	Description
Éteinte (sombre)	L'appareil est hors tension.
Lumière estompée	La porte de l'appareil est ouverte.
Clignotement lent jaune	L'alimentation de l'appareil est activée mais l'appareil n'est pas connecté au système de gestion de données chromatographiques.
Jaune	L'appareil est connecté au système de gestion de données chromatographiques.
Clignotement vert	L'appareil s'équilibre.
Verte	L'appareil est équilibré.
Bleue	Une injection ou une séquence est en cours d'exécution, y compris l'acquisition de données.
Rouge	Un problème ou une erreur est survenu. Pour obtenir le message correspondant, consultez l'Instrument Audit Trail. Pour connaître les mesures appropriées à prendre, reportez-vous à la section Dépannage (> page 133).

LED D'ÉTAT

La LED **STATUS** du panneau de commande, située à l'intérieur de l'appareil, fournit les informations suivantes :

LED D'ÉTAT	Description
Éteinte (sombre)	L'appareil est hors tension.
Verte	L'appareil fonctionne correctement.
Rouge	Un problème ou une erreur est survenu. Pour obtenir le message correspondant, consultez l'Instrument Audit Trail. Pour connaître les mesures appropriées à prendre, reportez-vous à la section Dépannage () page 133).

6.4 Mise sous et hors tension

L'interrupteur d'alimentation sur l'appareil est l'interrupteur principal permettant la mise sous et hors tension. L'interrupteur d'alimentation principal est placé en position Marche lors de l'installation initiale de l'appareil.

Afin de faciliter l'utilisation, vous pouvez utiliser le bouton d'alimentation situé sur la partie avant gauche du socle du système Vanquish (bouton d'alimentation du système) pour la mise sous et hors tension.

Respectez les consignes suivantes :

- Tous les modules du système Vanquish connectés au socle du système au moyen de câbles System Interlink sont mis sous ou hors tension en même temps lorsque vous appuyez sur le bouton d'alimentation du système.
- Lorsque le système est sous tension, le bouton d'alimentation du système est enfoncé. Lorsque le système est hors tension, le bouton d'alimentation du système est relâché.
- Si l'interrupteur d'alimentation principal d'un appareil est en position Arrêt, vous ne pouvez pas mettre cet appareil sous tension au moyen du bouton d'alimentation du système.
- Pour mettre un appareil complètement hors tension, vous devez placer son interrupteur d'alimentation principal en position Arrêt. Il ne suffit pas d'appuyer sur le bouton d'alimentation du système pour couper complètement l'alimentation électrique de l'appareil.

6.5 Préparation de l'appareil à l'utilisation

Cette section renseigne sur toutes les étapes supplémentaires requises pour préparer l'appareil à l'utilisation et à l'analyse d'échantillons.

Avant la toute première utilisation de l'appareil

Préparez l'appareil pour la première utilisation en prenant en compte de ce qui suit :

AVIS

Avant de commencer à utiliser l'appareil, rincez abondamment le trajet d'écoulement du système :

- Lorsque vous installez des appareils ou des composants dans le système, rincez-les et évacuez le liquide de rinçage avant de les raccorder au trajet d'écoulement du système. Pour rincer les modules Vanquish, suivez les instructions du Manuel d'utilisation du système Vanquish.
- Certains composants de l'appareil sont remplis d'isopropanol lors de l'expédition de celui-ci depuis le site de fabrication.Lorsque vous utilisez l'appareil pour la première fois, employez des solvants miscibles à l'isopropanol. À défaut, utilisez un solvant intermédiaire approprié.

Avant de commencer l'analyse d'un échantillon

Avant de débuter une analyse :

- Contrôlez le niveau de liquide dans les réservoirs de solvants. Vérifiez si la quantité de solvant est suffisante pour l'analyse.
- Fermez les portes de tous les modules du système Vanquish, si ce n'est pas déjà fait.
- Assurez-vous que le système chromatographique est correctement équilibré (voir plus bas).

Équilibrage du système

L'équilibrage du système doit comprendre les opérations suivantes :

- Purge de la pompe (*toutes* les voies, y compris celles non utilisées pour l'application)
- Rinçage de l'ensemble du système chromatographique avec le solvant initial pour éliminer tout solvant provenant d'une précédente analyse

- Mise à la température initiale (chauffage ou refroidissement) de tous les appareils du système équipés d'une régulation en température. Ce sont, par exemple :
 - le compartiment à colonnes et l'appareil de refroidissement post-colonne ;
 - le compartiment à échantillons thermostaté du le passeur d'échantillon ;
 - la cellule à écoulement d'un détecteur à fluorescence ;
 - le tube d'évaporation d'un détecteur d'aérosols chargés.
- Mise sous tension de la lampe (ou des lampes) dans un détecteur UV/VIS
- Surveillance de la pression de la pompe et de sa fluctuation ; vérification que la pression est stable et que la fluctuation est dans des limites raisonnables pour l'application
- Surveillance du signal du détecteur et contrôle de sa stabilité de telle sorte que la dérive et le bruit du signal restent dans des limites raisonnables pour l'application
- Exécution d'une mise à zéro automatique de la ligne de base du détecteur

NOTE Le logiciel Chromeleon prend en charge des procédures permettant de démarrer automatiquement un système chromatographique à partir du logiciel (**Smart Startup**). La procédure de démarrage couvre les opérations d'équilibrage du système. Pour obtenir plus de détails, reportez-vous à l'*Aide de Chromeleon*.

6.6 Principaux paramètres d'utilisation

Les paramètres décrits dans la présente section doivent être pris en considération pour l'utilisation routinière de l'appareil. Ces paramètres sont accessibles depuis l'interface utilisateur de Chromeleon.

Si l'un des paramètres répertoriés ci-dessous n'est pas disponible sous Chromeleon, pensez à mettre à jour la version du microprogramme et de Chromeleon.

Pour obtenir plus d'informations, reportez-vous à l'Aide et à la documentation Chromeleon.

NOTE Le Instrument Method Wizard (Assistant de méthode d'instrument) offre plusieurs vues différentes. Selon les paramètres à définir, vous pouvez choisir le mode de visualisation souhaité (par exemple **Easy** ou **Advanced**).

Paramètre	Description
ID de colonne	Le système d'identification des colonnes (ID de colonne) vous permet de stocker des informations spécifiques sur un connecteur d'ID de colonne. Pour obtenir une liste complète des propriétés des colonnes et des renseignements complémentaires sur le système d'identification de colonne, reportez-vous à la section Identification des colonnes (ID de colonne) (> page 35) ou consultez l'Aide de Chromeleon. Pour activer l'identification de colonne sous Chromeleon,
	reportez-vous à la section Installation de l'identification des colonnes (ID de colonne) (▶ page 66).
Sélection de colonne	Si le compartiment à colonnes est doté d'une ou de deux vannes de commutation pour colonne, la sélection de colonne s'effectue sous Chromeleon. Le nombre de vannes de commutation pour colonne, le modèle de vanne et l'application déterminent les commandes et les propriétés à employer.
Equilibration Time	Le paramètre Equilibration Time peut être défini pour le compartiment à colonnes, l'appareil de refroidissement post-colonne (VH-C10 uniquement) et les appareils de préchauffage actifs (s'ils sont installés).
	Pour obtenir plus de détails, reportez-vous à la section Delta de température de préparation et durée d'équilibrage () page 93) ou consultez l'Aide de Chromeleon.

Paramètre	Description
Fast Cool	Par défaut, le mode Fast Cool est désactivé. Pour refroidir rapidement les surfaces chaudes situées à l'intérieur du compartiment à colonnes, activez le mode Fast Cool sous Chromeleon ou appuyez sur le bouton FAST COOL du panneau de commande. Reportez-vous à la section Panneau de commande () page 84).
	Pour annuler le mode Fast Cool sous Chromeleon :
	 Definissez une nouvelle temperature nominale pour l'appareil de refroidissement post-colonne, l'appareil de préchauffage actif ou le compartiment à colonnes. Le composant concerné est chauffé jusqu'à atteindre la température de consigne. Notez que la régulation de la température reste désactivée pour les composants pour lesquels aucune nouvelle température nominale n'a été définie.
	 Activez la régulation de la température pour l'appareil de refroidissement post-colonne, l'appareil de préchauffage actif ou le compartiment à colonnes en cliquant sur ON. Notez que la régulation en température reste désactivée pour les composants pour lesquels la régulation de température est désactivée.
Paramètres des capteurs de fuites	Par défaut, le capteur de fuites est activé.
Ready Temp Delta	Le paramètre Ready Temp Delta peut être réglé pour le compartiment à colonnes (VH-C10 uniquement), l'appareil de refroidissement post-colonne et les appareils de préchauffage actifs (s'ils sont installés).
	Pour obtenir plus de détails, reportez-vous à la section Delta de température de préparation et durée d'équilibrage (▶ page 93) ou consultez l'Aide de Chromeleon.
Thermostatisation des colonnes	Dans la fenêtre Command , sélectionnez le compartiment à colonnes, nommé par défaut CC , puis sélectionnez ON pour TempCtrl pour activer la thermostatisation des colonnes. Pour modifier la température de thermostatisation, saisissez une valeur pour le paramètre Temperature.Nominal .
Thermostatisation de l'appareil de refroidissement post- colonne (VH-C10 uniquement)	Dans la fenêtre Command , sélectionnez l'appareil de refroidissement post-colonne, nommé par défaut PCC , puis sélectionnez ON pour TempCtrl pour activer la thermostatisation de l'appareil de refroidissement post- colonne.
	Pour modifier la température de thermostatisation pour l'appareil de refroidissement post-colonne, saisissez une valeur pour le paramètre Temperature.Nominal .

Paramètre	Description
Thermostatisation des appareils de préchauffage actifs (s'ils sont installés)	Dans la fenêtre Command , sélectionnez les appareils de préchauffage actifs, nommé par défaut PrehLeft et PrehRight , puis sélectionnez ON pour TempCtrl pour activer la thermostatisation des appareils de préchauffage actifs. Pour modifier la température de thermostatisation pour les appareils de préchauffage actifs, saisissez une valeur pour le paramètre Temperature.Nominal .
Modes de thermostatisation	Le mode air calme est sélectionné par défaut. Si nécessaire, modifiez le paramètre et sélectionnez le mode Forced air. Pour obtenir des renseignements complémentaires sur les modes de thermostatisation, reportez-vous à la section Principe de fonctionnement () page 31).
Signaux de température	Dans la boîte de dialogue Instrument Configuration Manager pour le compartiment à colonnes, les cases de température sont cochées par défaut lorsque le compartiment à colonnes est configuré sous Chromeleon. Le logiciel Chromeleon utilise ce paramétrage pour générer les canaux d'enregistrement des signaux de température.
	Si un problème survient, les signaux de température peuvent fournir des renseignements utiles à l'identification et à l'élimination de la source du problème. Par conséquent, enregistrez toujours les signaux de température.

6.7 Delta de température de préparation et durée d'équilibrage

La propriété **Ready** indique si le compartiment à colonnes est prêt à fonctionner. Lorsque l'état du compartiment à colonnes est **NotReady**, vous ne pouvez pas lancer l'analyse. Gardez à l'esprit que les paramètres **Ready Temp Delta** et **Equilibration Time** influencent la durée nécessaire au compartiment à colonnes pour passer à l'état **Ready**.

Le logiciel Chromeleon prédéfinit le paramètre **Equilibration Time** par défaut. Ce paramètre peut être ajusté si besoin.

Ready Temp Delta

Le paramètre **Ready Temp Delta** indique l'écart positif ou négatif toléré par rapport à la température de consigne. Si l'écart entre la température actuelle et la valeur de consigne est supérieur à la valeur saisie ici, le compartiment à colonnes passe à l'état **NotReady** et n'est pas prêt à fonctionner.

Exemple : température de consigne = 45°C, Ready Temp Delta = 1°C

Le compartiment à colonnes est prêt à fonctionner si la température actuelle du paramètre **Equilibration Time** est comprise entre 44 et 46°C. Si la température actuelle est en dehors de cette plage, l'état du compartiment à colonnes est **NotReady**.

Temps d'équilibrage

Ce paramètre indique la durée pendant laquelle la température souhaitée doit être dans la plage définie par la température de consigne et le paramètre de **Ready Temp Delta** avant que le compartiment à colonnes ne soit prêt pour le fonctionnement et que l'analyse ne puisse démarrer.

Exemple : température de consigne = 45°C, **Ready Temp Delta** = 1°C, **Equilibration Time** = 0,5 min

Le compartiment à colonnes est prêt à fonctionner si la température actuelle est comprise entre 44 et 46°C et si elle est restée dans cette plage pendant 0,5 minute.

Respectez les consignes suivantes :

 Plus la valeur de Ready Temp Delta est faible et plus la valeur de Equilibration Time est élevée, plus la durée nécessaire avant que le compartiment à colonnes ne soit prêt à fonctionner et avant que l'acquisition des données et/ou l'analyse ne puissent démarrer est importante.

- Les conditions de l'environnement de travail, par exemple un courant d'air ou une climatisation, peuvent également augmenter la durée d'attente voire empêcher que le compartiment à colonnes ne passe à l'état **Ready**.
- Cependant, gardez à l'esprit que plus le système est bien équilibré, meilleurs sont les résultats de chromatographie.

6.8 Optimisation des performances de l'appareil

Cette section fournit des informations sur les performances optimales de l'appareil et vous indique comment améliorer encore les performances.

6.8.1 Consignes générales

Pour optimiser les performances du compartiment à colonnes, suivez les consignes générales ci-après :

- Utilisez des capillaires de raccordement courts.
- Surveillez l'utilisation des composants spécifiques du compartiment à colonnes, par exemple les vannes de commutation pour colonne sujettes à l'usure, et prévoyez une périodicité adéquate des opérations de maintenance (reportez-vous à la section Fonctions Predictive Performance (> page 111)).
- Fermez toujours le capot de chambre à colonnes et la porte du compartiment à colonnes pendant le fonctionnement.
- Respectez les instructions et recommandations générales sur l'utilisation de solvants et d'additifs dans le système chromatographique. Consultez la section Utilisation de solvants et additifs du Manuel d'utilisation du système Vanquish.

6.8.2 Préchauffage actif

Un appareil de préchauffage actif peut être installé afin de porter l'éluant à la température souhaitée avant l'analyse. Pour obtenir des informations concernant l'installation d'un appareil de préchauffage actif, reportez-vous à la section Raccordement et installation de l'appareil de préchauffage actif (> page 73).

6.9 Arrêt de l'appareil

Si une période d'inactivité de l'appareil est prévue, suivez les instructions de cette section pour le mettre à l'arrêt.

NOTE Le logiciel Chromeleon fournit des procédures automatiques de préparation à l'arrêt du système de chromatographie. Ces procédures couvrent notamment les opérations de diminution du débit, de réduction de la température des appareils à température régulée et de mise hors tension des lampes du détecteur. Pour plus d'informations sur les fonctionnalités **Smart Shutdown** et **Smart Standby**, consultez l'*Aide de Chromeleon*.

6.9.1 Arrêt de courte durée (interruption de fonctionnement)

Pour interrompre le fonctionnement de l'appareil pendant une courte durée (arrêt de courte durée), par exemple la nuit, suivez les instructions suivantes, en fonction des modules présents dans votre système Vanquish :

• Concernant votre détecteur Vanquish, veuillez noter les points suivants :

Type de détecteur	Description
Détecteur d'aérosols chargés	Vérifiez qu'il y a suffisamment de gaz pour conserver un débit de gaz à travers le détecteur. Ceci permet d'empêcher toute formation de résidus provenant des solvants ou analytes. Le gaz doit circuler lors du refoulement du débit de la pompe vers le détecteur.
Détecteurs UV/VIS	La ou les lampes du détecteur peuvent rester allumées. <i>Pour les détecteurs à longueur d'onde variable et à barrette de diodes VH-D10 seulement :</i> L'obturateur peut être placé en position fermée pour protéger la cellule à écoulement.
Détecteur à fluorescence	Coupez la régulation de température pour la cellule à écoulement.

- Définissez un débit de 0,05 mL/min, puis réglez la pompe de sorte que le solvant refoulé soit approprié.
 Vérifiez la limite de pression inférieure de la pompe, et modifiez la valeur si nécessaire. Si la pression chute en dessous de la limite inférieure, la pompe interrompt le débit.
- Mettez la vanne d'injection du passeur d'échantillon en position d'injection (Inject).
- Assurez-vous que la température de la colonne ne dépasse pas 40 °C.

 Lors de la reprise du fonctionnement, laissez le débit s'équilibrer puis, avant de continuer, vérifiez que les paramètres de fonctionnement des autres modules du système sont définis conformément à vos besoins.

6.9.2 Arrêt de longue durée

Pour interrompre le fonctionnement pour une période prolongée, suivez les instructions ci-dessous.

NOTE L'arrêt de l'appareil a un effet sur le fonctionnement du système. Lors de l'arrêt de l'appareil, respectez également les instructions de mise à l'arrêt des autres modules du système Vanquish et appliquez les mesures appropriées (consultez les *Manuels d'utilisation* des autres modules concernés).

- 1. Retirez la colonne.
- 2. Rincez le système avec un solvant adéquat (au minimum de qualité HPLC). Respectez les consignes suivantes :

NOTE Lorsqu'un système Vanquish Core a été modifié pour utiliser des additifs et des solvants compatibles avec la phase normale, consultez les informations concernant le liquide de rinçage dans la section *Considérations relatives aux additifs et solvants compatibles avec la phase normale* du *Manuel d'utilisation* du *système Vanquish*.

Situation après arrêt	Si aucun additif n'est utilisé	Si un additif a été utilisé
L'appareil reste dans le laboratoire après l'arrêt	Rincez le système avec du méthanol, par exemple. N'utilisez pas de l'acétonitrile à 100 %.	Rincez le système avec plusieurs volumes de méthanol et d'eau (50:50) (par exemple, 1,0 mL/min pendant 10 minutes pour le système standard) pour éviter toute accumulation de sels dans le circuit fluidique. Si les solvants employés dans l'appareil ne sont pas miscibles à l'eau, utilisez un solvant intermédiaire approprié.

Situation après arrêt	Si aucun additif n'est utilisé	Si un additif a été utilisé
L'appareil doit être transporté ou expédié après l'arrêt	Rincez le système avec de l'isopropanol.	Rincez d'abord le système avec plusieurs volumes de méthanol et d'eau (50:50) (par exemple, 1,0 mL/min pendant 10 minutes pour le système standard) pour éviter toute accumulation de sels dans le circuit fluidique. Si les solvants employés dans l'appareil ne sont pas miscibles à l'eau, utilisez un solvant intermédiaire approprié. Ensuite, rincez le système avec de l'isopropanol.

- 3. Arrêtez le débit de la pompe. Patientez jusqu'à ce que la pression du système redescende à zéro avant de poursuivre l'arrêt de l'appareil.
- 4. Cette étape dépend des critères ci-après :

Situation	Étapes
L'appareil et tous les modules du système restent dans le système empilé et vont être arrêtés	Arrêtez le système au moyen du bouton d'alimentation électrique situé sur le socle du système.
L'appareil doit être transporté ou expédié après l'arrêt	Si l'un des modules doit être retiré du système empilé, arrêtez <i>tous</i> les modules du système à l'aide de leur interrupteur d'alimentation principal. Il ne suffit pas d'appuyer sur le bouton d'alimentation du système pour couper complètement l'alimentation électrique des appareils.
	Suivez les instructions sous Déménagement ou expédition de l'appareil () page 129).

6.9.3 Redémarrage après un arrêt de longue durée

Pour redémarrer l'appareil après un arrêt de longue durée, suivez les étpaes suivantes :

1. Mettez l'appareil sous tension. Respectez les consignes suivantes :

Situation	Action
Si l'appareil est resté dans le système empilé et tous les modules du système sont hors tension.	Mettez le système sous tension au moyen du bouton d'alimentation électrique situé sur le socle du système.
Si l'appareil est redémarré après le transport.	Mettez l'appareil sous tension au moyen de l'interrupteur d'alimentation principal.

- 2. Préparez et redémarrez les autres modules du système Vanquish en suivant les instructions des *Manuels d'utilisation* de chaque module. Veuillez accorder une attention particulière à la section *Préparation au fonctionnement du module*.
- 3. Avant de lancer une analyse, laissez l'appareil s'équilibrer et assurezvous qu'il est prêt à fonctionner. Consultez la section Préparation de l'appareil à l'utilisation (▶ page 88).

7 Maintenance et entretien

Ce chapitre décrit les procédures de maintenance et d'entretien de routine que l'utilisateur peut être amené à effectuer.

7.1 Présentation de la maintenance et de l'entretien

Ce chapitre décrit les procédures de maintenance, d'entretien et de réparation de routine que l'utilisateur peut être amené à effectuer.



Seul le personnel d'entretien certifié par Thermo Fisher Scientific (par souci de concision, désigné par la suite par techniciens d'entretien Thermo Fisher Scientific) est autorisé à réaliser des procédures supplémentaires de maintenance et d'entretien.

L'appareil est conçu de manière à faciliter la maintenance et l'entretien. Les parties de l'appareil pouvant être entretenues par l'utilisateur sont accessibles depuis l'avant. Sauf mention contraire, les procédures de maintenance ne nécessitent pas de retirer l'appareil du système.

Les procédures de maintenance ne nécessitent pas de retirer la porte. Toutefois, si cela est nécessaire pour quelque raison que ce soit ou dans le cadre d'une procédure spécifique, il est possible de retirer une porte. Si vous devez retirer une porte, suivez les étapes figurant dans la section Remplacement de la porte de l'appareil () page 125).

7.2 Consignes de sécurité pour la maintenance et l'entretien

Lorsque vous effectuez des procédures de maintenance ou d'entretien, respectez les consignes de sécurité suivantes :



Respectez tous les messages d'avertissement et mentions de mise en garde figurant à la section Précautions de sécurité (> page 19).



AVERTISSEMENT—Haute tension

Des tensions élevées sont présentes dans l'instrument et peuvent provoquer des décharges électriques.

N'ouvrez pas le boîtier et ne retirez pas les capots de protection, à moins que cela ne soit expressément indiqué dans le présent manuel.



AVERTISSEMENT—Fuite de substances dangereuses provenant des raccordements fluidiques

Les raccordements fluidiques et les connexions capillaires peuvent contenir des substances potentiellement dangereuses pour la santé. Il existe un risque de projection de solvant si des capillaires éclatent, se libèrent de leurs raccords, ou si les raccords ne sont pas suffisamment serrés ou se déconnectent.

- Portez un équipement de protection approprié et suivez les bonnes pratiques de laboratoire.
- Avant d'entamer les procédures de maintenance ou de réparation, évacuez toutes les substances nocives à l'aide d'un solvant approprié.



AVERTISSEMENT—Inclinaison des réservoirs de liquides

Les liquides contenus dans les réservoirs placés dans le bac à solvants peuvent contenir des substances nocives. Ces substances peuvent présenter des risques pour la santé et la sécurité si elles se répandent.

Afin d'éviter tout déversement provenant des réservoirs, faites attention de ne pas tirer sur les conduites des liquides lors de la maintenance.



ATTENTION—Projection de solvant

Un risque de projection de solvants existe lorsque ceux-ci sont soumis à une pression élevée.

- Arrêtez le débit de la pompe avant ouverture du trajet d'écoulement.
- Patientez jusqu'à ce que la pression du système redescende à zéro.
- Revêtez un équipement de protection approprié lorsque vous ouvrez le trajet d'écoulement.



ATTENTION—Surfaces chaudes

Les surfaces internes de l'appareil peuvent s'échauffer lors de l'opération. Ne touchez pas les surfaces chaudes pour éviter de vous brûler.

- Ne touchez pas la face interne du capot de la chambre à colonnes. Lorsque vous ouvrez le capot de la chambre à colonnes, saisissez-le uniquement par le porte-capot.
- Attendez que les surfaces chaudes aient refroidi avant de commencer toute procédure de remplacement ou de maintenance.



ATTENTION—Pression hydrostatique

Il existe un risque de déversement de solvants lorsque vous ouvrez le trajet d'écoulement. Cela est dû à la pression hydrostatique présente dans le système lorsque les réservoirs de solvants sont situés au-dessus de l'orifice de refoulement de la pompe. Avant de desserrer un raccordement dans le trajet d'écoulement :

- Arrêtez le débit de la pompe et patientez jusqu'à ce que la pression du système redescende à zéro.
- Dévissez les bouchons des réservoirs de solvants et retirez les conduites de solvant en même temps que les bouchons des réservoirs.
- Videz les conduites de solvant. Pour obtenir plus de détails, consultez le Manuel d'utilisation de la pompe.
- Ensuite, resserrez les bouchons des réservoirs.



ATTENTION—Décharge électrique ou détérioration de l'appareil

Lorsque l'appareil est éteint, il n'est pas complètement isolé électriquement tant que le cordon d'alimentation est branché. Réaliser des réparations alors que l'appareil est connecté à l'alimentation électrique peut entraîner des accidents corporels.

- Débranchez toujours le cordon d'alimentation avant de procéder à des travaux de réparation à l'intérieur de l'appareil.
- Si vous devez enlever des capots ou des panneaux, ne branchez pas le cordon d'alimentation à l'appareil tant que les panneaux et les capots sont démontés.

7.3 Règles générales de maintenance et d'entretien

Conformez-vous aux règles et aux recommandations suivantes afin de procéder correctement à la maintenance et à l'entretien.

Règles générales

- Avant de débuter la procédure de maintenance ou d'entretien, arrêtez l'appareil lorsque vous en recevez l'instruction.
- Employez exclusivement les pièces de rechange spécifiquement autorisées et certifiées pour l'appareil par Thermo Fisher Scientific.
- Suivez toutes les instructions étape par étape et utilisez les outils recommandés pour la procédure.

Utilisation du mode Fast Cool pour les procédures de maintenance

Pour refroidir rapidement les surfaces chaudes situées à l'intérieur du compartiment à colonnes, utilisez le mode Fast Cool :

- Appuyez sur le bouton FAST COOL du panneau de commande. Lorsque la LED à côté du bouton devient verte, vous pouvez accéder au compartiment à colonnes.
- Annulez le mode Fast Cool une fois la procédure terminée, en appuyant de nouveau sur le bouton FAST COOL. Annulez le mode Fast Cool une fois la procédure terminée via le logiciel Chromeleon (reportez-vous à la section Principaux paramètres d'utilisation (> page 90)).

Ouverture des connexions du trajet d'écoulement

- Avant d'ouvrir le trajet d'écoulement afin de remplacer les capillaires du système, arrêtez la pompe et patientez jusqu'à ce que la pression du système redescende à zéro.
- Les composants sales peuvent contaminer le système chromatographique. La contamination diminue les performances des modules et du système dans son ensemble, ou peut même causer une détérioration des modules et du système. Par conséquent :
 - Portez toujours des gants appropriés.
 - Placez les composants uniquement sur une surface propre et non pelucheuse.
 - Veillez à maintenir vos outils propres.
 - Utilisez uniquement des chiffons non pelucheux pour le nettoyage.

Réparations en atelier

 Si vous devez renvoyer l'appareil pour une réparation au dépôt, suivez les instructions figurant dans la section Déménagement ou expédition de l'appareil (> page 129).

7.4 Maintenance de routine et maintenance préventive

Seul un appareil en bon état et correctement entretenu vous permettra d'optimiser les performances, la disponibilité de l'appareil et l'exactitude des résultats.

7.4.1 Plan de maintenance

Effectuez régulièrement les procédures de maintenance figurant dans le tableau ci-dessous. Les fréquences indiquées dans le tableau ne sont que des suggestions. La fréquence optimale des opérations de maintenance dépend de plusieurs facteurs, tels que le type et la quantité d'échantillons et de solvants utilisés avec l'appareil.

Fréquence	Opération de maintenance à réaliser
Tous les jours	 Inspectez les raccordements fluidiques afin d'y détecter d'éventuels signes de fuite ou d'obstruction. Serrez les raccords qui présentent une fuite. Si vous utilisez des tampons ou des solutions salines, rincez soigneusement l'appareil après utilisation au moyen d'un solvant adéquat ne contenant ni tampon ni sel.
Régulièrement	 Inspectez les raccordements fluidiques afin d'y détecter d'éventuels dommages, comme des fissures, des entailles, des coupures ou des obstructions. Le tube d'évacuation est raccordé à l'orifice d'évacuation situé dans la partie avant basse de l'appareil, au-dessous de la porte avant. Vérifiez que le tube n'est pas bouché et qu'il est bien acheminé sous l'orifice d'évacuation. Si nécessaire, videz le bac d'évacuation. Vérifiez que toutes les étiquettes d'avertissement sont bien présentes sur l'appareil et qu'elles sont lisibles. Dans le cas contraire, prenez contact avec Thermo Fisher Scientific pour un remplacement.
Annuellement	Demandez au personnel d'entretien de Thermo Fisher Scientific de réaliser une maintenance préventive une fois par an.

NOTE Le logiciel Chromeleon est doté de fonctions permettant d'estimer la durée de vie des consommables (reportez-vous à la section Fonctions Predictive Performance () page 111).
7.4.2 Nettoyage ou décontamination de l'appareil

Le nettoyage et la décontamination doivent être réalisés par du personnel qualifié portant des équipements de protection individuelle appropriés. Respectez toujours les réglementations nationales et locales.

AVIS

Essuyez immédiatement tout liquide renversé sur le système. Une exposition prolongée à ces liquides pourrait endommager les surfaces.

Décontamination

Une décontamination est par exemple requise en présence d'une fuite ou d'un déversement accidentel, ou avant de procéder à l'entretien ou au transport de l'appareil. Utilisez un détergent ou désinfectant approprié afin de vous assurer que le traitement permette une manipulation sans danger de l'appareil.

Éléments nécessaires

- Détergent approprié (ou désinfectant)
- Eau purifiée
- Chiffon ou serviettes en papier non pelucheux



ATTENTION—Mélanges de gaz explosifs provenant de détergents à base d'alcool

Les détergents contenant de l'alcool peuvent former des mélanges de gaz inflammables et explosifs s'ils sont exposés à l'air.

- Utilisez ces détergents uniquement dans les cas requis et uniquement dans des pièces bien ventilées.
- Évitez la proximité des flammes nues ou toute exposition à une chaleur excessive durant le processus de nettoyage.
- Essuyez les composants nettoyés jusqu'à ce qu'ils soient secs après nettoyage. N'utilisez pas l'appareil tant qu'elle n'est pas entièrement sèche.

AVIS

Respectez les consignes suivantes :

- Utilisez uniquement des détergents qui n'endommagent pas les surfaces du système.
- N'utilisez jamais d'outils tranchants ou de brosses pour nettoyer les surfaces.
- N'utilisez pas de vaporisateurs pour le nettoyage.
- Évitez que du détergent ne pénètre dans le trajet d'écoulement.
- Le chiffon ou les serviettes en papier employés pour le nettoyage ne doivent pas être trop humides. Évitez que du liquide ne pénètre dans les composants fonctionnels de l'appareil. Les liquides peuvent provoquer un court-circuit en entrant en contact avec les composants électroniques.

Préparatifs

1. Mettez l'appareil hors tension, puis débranchez le cordon d'alimentation de la source d'alimentation.

- Essuyez les surfaces au moyen d'un chiffon ou d'une serviette en papier propre, sec, doux et non pelucheux. Si nécessaire, humectez légèrement le chiffon ou la serviette en papier avec une solution d'eau tiède et de détergent approprié.
- 2. Laissez le détergent agir conformément aux recommandations du fabricant.
- 3. Essuyez les surfaces nettoyées avec de l'eau purifiée afin d'éliminer toute trace de détergent.
- 4. Séchez les surfaces au moyen d'un chiffon ou d'une serviette en papier doux et non pelucheux.

7.4.3 Fonctions Predictive Performance

Le logiciel Chromeleon est doté de fonctions permettant de suivre et d'enregistrer les informations d'entretien et de qualification de l'appareil.

Surveillance des intervalles d'entretien et de qualification

Les panneaux dédiés à l'entretien et à la qualification vous permettent de définir la périodicité des procédures d'entretien ou de qualification. Ces fonctions, appelées Predictive Performance, vous permettent de programmer ces procédures sur la base des conditions réelles de fonctionnement et d'utilisation de l'appareil. Vous pouvez également définir des limites afin d'être prévenu de la date d'entretien ou de qualification.

Les barres à code-couleurs offrent un retour visuel qui vous permet de consulter et de surveiller facilement l'état de l'appareil. Si une limite d'avertissement a été définie, un message affiché dans la Chromeleon Audit Trail vous informe lorsqu'une action est prévue.

Les compteurs d'entretien et de qualification peuvent être remis à zéro une fois l'action requise effectuée. Pour maintenir à jour les informations de Predictive Performance, veillez à remettre à zéro le compteur dès qu'une procédure d'entretien ou de qualification a été effectuée.

Pour obtenir plus d'informations, reportez-vous à l'Aide de Chromeleon.

7.5 Vanne de commutation pour colonne

Le compartiment à colonnes peut accueillir une ou deux vannes de commutation pour colonne. Pour assurer un fonctionnement optimal de la vanne de commutation pour colonne, suivez les consignes ci-après :

- L'utilisation, l'entreposage ou le transport de la vanne à une température inférieure à 0°C avec de l'eau dans les trajets d'écoulement peut endommager les surfaces d'étanchéité.
- Installez des raccords du connecteur Viper pour le raccordement des capillaires à la vanne.
- Nettoyez la vanne en rinçant toutes les conduites à l'aide d'un solvant adéquat. La nature du solvant à employer dépend des échantillons et des phases mobiles utilisées.

7.5.1 Inspection de la vanne de commutation pour colonne à la recherche de fuites

À quel moment exécuter cette opération ?

- Des gouttelettes de liquide apparaissent au niveau de l'orifice de la vanne
- Du liquide s'est accumulé dans le bac de fuite
- Des gouttelettes de liquide apparaissent autour de l'orifice d'évacuation



Illustration 35: L'orifice d'évacuation sur la vanne de commutation pour colonne

Procédez comme suit

1. Comme les fuites surviennent généralement aux raccords, inspectez visuellement tous les raccords du trajet d'écoulement.

Situation	Action
Des gouttelettes de liquide apparaissent au niveau de l'orifice de la vanne	 Resserez le capillaire affecté. Si des gouttelettes apparaissent toujours au niveau de l'orifice de la vanne, changez le capillaire.
Du liquide s'est accumulé dans le bac de fuite	 Reserrez tous les capillaires. Si le liquide continue de s'accumuler dans le bac de fuite, changez les capillaires.
Des gouttelettes de liquide apparaissent autour de l'orifice d'évacuation	 Si c'est le cas, absorbez-les avec une serviette en papier. Si des gouttelettes continuent à apparaître autour de l'orifice d'évacuation, remplacez la vanne de commutation pour colonne. Reportez-vous à la section Remplacement de la vanne de commutation pour colonne (> page 113).

2. Cette étape dépend des critères ci-après :

7.5.2 Remplacement de la vanne de commutation pour colonne

À quel moment exécuter cette opération ?

- La vanne de commutation pour colonne fuit
- Pour changer entre différents types de vanne

Éléments nécessaires

Vanne de commutation pour colonne, selon les besoins de l'application

Outils nécessaires

Tournevis type Torx T10

Préparatifs



AVERTISSEMENT—Fuite de substances dangereuses provenant des raccordements fluidiques

Les raccordements fluidiques et les connexions capillaires peuvent contenir des substances potentiellement dangereuses pour la santé. Il existe un risque de projection de solvant si des capillaires éclatent, se libèrent de leurs raccords, ou si les raccords ne sont pas suffisamment serrés ou se déconnectent.

- Portez un équipement de protection approprié et suivez les bonnes pratiques de laboratoire.
- Avant d'entamer les procédures de maintenance ou de réparation, évacuez toutes les substances nocives à l'aide d'un solvant approprié.

- 1. Arrêtez le débit de la pompe.
- 2. Débranchez toutes les conduites de liquide raccordées à la vanne de commutation pour colonne.
- 3. Désactivez la vanne dans la boîte de dialogue Instrument Configuration Manager.
- Mettez le compartiment à colonnes hors tension. S'il est sous tension, la vanne de commutation peut être endommagée pendant le remplacement.
- Ne retirez pas encore les bouchons de raccordement qui couvrent les orifices des vannes sur la vanne de commutation pour colonne de remplacement.

Retrait de la vanne de commutation pour colonne

- 1. Desserrez les 4 vis de la vanne de commutation pour colonne.
- 2. Conservez les vis pour la pose de la vanne de commutation pour colonne de remplacement.



Illustration 36: Retrait de la vanne de commutation pour colonne

- 3. Retirez avec précaution du boîtier la vanne de commutation pour colonne.
- 4. Déconnectez le câble de l'actionneur de la vanne situé à l'arrière de la vanne de commutation pour colonne.
- S'il n'est pas prévu d'installer ultérieurement une vanne de commutation, remettez en place le capot par dessus l'emplacement pour la vanne de commutation pour colonne. Reportez-vous à la section Remplacement du capot de la vanne de commutation pour colonne (▶ page 116).

Installation de la vanne de commutation pour colonne

- 1. Contrôlez l'orientation du câble. La broche d'alignement située au milieu du connecteur du câble doit être dirigée vers le bas.
- 2. Branchez le câble au connecteur situé à l'arrière de la vanne de commutation pour colonne.
- 3. Enfoncez le câble dans le boîtier.
- 4. Enfoncez délicatement la vanne de commutation pour colonne dans le boîtier.





- 5. À l'aide du tournevis, serrez les 4 vis de la vanne de commutation pour colonne.
- 6. Retirez les bouchons de raccordement qui couvrent les orifices des vannes.
- Connectez les conduites de liquide, selon les besoins de l'application. Pour les instructions, reportez-vous à la section Raccordement des composants à une vanne de commutation pour colonne (▶ page 76).
- 8. Mettez le compartiment à colonnes sous tension.
- 9. Configurez la vanne de commutation pour colonne dans le système de gestion de données chromatographiques :
 - a) Dans la boîte de dialogue Instrument Configuration Manager, sélectionnez l'emplacement (**Upper Valve**, **Lower Valve**).
 - b) Sélectionnez le modèle de vanne.
- 10. Mettez la vanne dans les différentes positions, puis testez les raccordements fluidiques.

7.5.3 Remplacement du capot de la vanne de commutation pour colonne

Éléments nécessaires

Capot de vanne de commutation pour colonne

Outils nécessaires

Tournevis type Torx T10

Retrait du capot de la vanne de commutation pour colonne

- 1. Ouvrez la porte du compartiment à colonnes.
- 2. À l'aide du tournevis, desserrez les 4 vis du capot de la vanne de commutation.



Illustration 38: Vis du capot de la vanne de commutation

- 3. Retirez les vis et conservez-les pour une installation ultérieure.
- 4. Retirez le capot et le joint d'étanchéité de la vanne de commutation.



Illustration 39: Retrait du capot et du joint de la vanne de commutation

N°	Description
1	Joint
2	Capot

5. Détachez avec précaution le câble de l'arrière du joint d'étanchéité pour vanne de commutation.

Installation du capot de vanne de commutation pour colonne

- 1. Fixez le câble à l'arrière du nouveau capot/joint d'étanchéité pour vanne de commutation pour colonne.
- 2. Placez le joint d'étanchéité et le capot par-dessus l'emplacement pour la vanne de commutation pour colonne.
- 3. À l'aide du tournevis, serrez les 4 vis sur le capot de vanne de commutation pour colonne.

7.6 Remplacement de l'appareil de refroidissement post-colonne (VH-C10 uniquement)

À quel moment exécuter cette opération ?

L'appareil de refroidissement post-colonne est défectueux

Éléments nécessaires

- Appareil de refroidissement post-colonne
- 2 vis

Préparatifs



ATTENTION—Surfaces chaudes

Les surfaces internes de l'appareil peuvent s'échauffer lors de l'opération. Ne touchez pas les surfaces chaudes pour éviter de vous brûler.

- Laissez-le refroidir avant de commencer l'opération d'entretien.
- Ne touchez pas la face interne du capot de la chambre à colonnes. Lorsque vous ouvrez le capot de la chambre à colonnes, saisissez-le uniquement par le porte-capot.
- Attendez que les surfaces chaudes aient refroidi avant de commencer toute procédure de remplacement ou de maintenance.

NOTE Pour refroidir rapidement l'appareil de refroidissement postcolonne, appuyez sur le bouton **FAST COOL** du panneau de commande. Patientez jusqu'à ce que la LED du bouton soit verte.

- 1. Ouvrez la porte du compartiment à colonnes.
- Débranchez les capillaires de l'appareil de refroidissement postcolonne du détecteur et de la vanne de commutation pour colonne, ou de la colonne.

3. Desserrez les 2 vis de l'appareil de refroidissement post-colonne.



Illustration 40: Retrait des vis de l'appareil de refroidissement postcolonne

- 4. Retirez l'appareil de refroidissement post-colonne de la plaque de montage.
- Pour monter le nouvel appareil de refroidissement post-colonne, suivez les instructions de la section Installation de l'appareil de refroidissement post-colonne (VH-C10 uniquement) (▶ page 59).

NOTE Annulez le mode Fast Cool une fois la procédure terminée, en appuyant de nouveau sur le bouton **FAST COOL**.

7.7 Remplacement de l'appareil de préchauffage

Cette section traite du remplacement d'un appareil de préchauffage dans le compartiment à colonnes. Selon la configuration du compartiment à colonnes, différents types d'appareils de préchauffage sont disponibles.

- Appareil de préchauffage actif, reportez-vous à la section Remplacement de l'appareil de préchauffage actif (> page 120).
- Appareil de préchauffage passif, reportez-vous à la section Remplacement de l'appareil de préchauffage passif (> page 122).

7.7.1 Remplacement de l'appareil de préchauffage actif

À quel moment exécuter cette opération ?

L'appareil de préchauffage actif est défectueux

Éléments nécessaires

Appareil de préchauffage actif

Préparatifs



ATTENTION—Surfaces chaudes

Le compartiment à colonnes et l'appareil de préchauffage peuvent être chauds. Ne touchez pas les surfaces chaudes pour éviter de vous brûler. Laissez-les refroidir avant de commencer l'opération d'entretien.

- 1. Détachez le porte-colonne du raccord de l'appareil de préchauffage actif.
- 2. Débranchez l'appareil de préchauffage actif de la colonne.



Illustration 41: Démontage de l'appareil de préchauffage actif de la colonne

3. Débranchez le câble de l'appareil de préchauffage du port situé à côté de la vanne de commutation inférieure.



Illustration 42: Port de connexion de l'appareil de préchauffage

- 4. Retirez le câble de l'appareil de préchauffage du serre-câble situé au-dessus du port de préchauffage.
- 5. Remplacez l'appareil de préchauffage défectueux.
- Raccordez le nouvel appareil de préchauffage à la colonne. Pour les instructions, reportez-vous à la section Raccordement et installation de l'appareil de préchauffage actif (▶ page 73).

7.7.2 Remplacement de l'appareil de préchauffage passif

À quel moment exécuter cette opération ?

L'appareil de préchauffage est défectueux

Éléments nécessaires

Appareil de préchauffage passif

Outils nécessaires

Tournevis type Torx T10

Procédez comme suit



ATTENTION—Surfaces chaudes

Le compartiment à colonnes et l'appareil de préchauffage peuvent être chauds. Ne touchez pas les surfaces chaudes pour éviter de vous brûler. Laissez-les refroidir avant de commencer l'opération d'entretien.

- 1. Débranchez le capillaire de sortie de l'appareil de préchauffage passif de l'entrée de la colonne.
- Débranchez le capillaire d'entrée de l'appareil de préchauffage passif du passeur d'échantillon ou de la vanne de commutation pour colonne.
- 3. À l'aide du tournevis, desserrez les deux vis et rondelles sur l'appareil de préchauffage passif et retirez-le.
- Installez le nouvel appareil de préchauffage dans la chambre à colonnes. Reportez-vous à la section Installation de l'appareil de préchauffage passif (▶ page 61).

7.8 Remplacement des fusibles de l'alimentation principale

À quel moment exécuter cette opération ?

Fusibles grillés

Éléments nécessaires

Fusibles (2 fusibles, 4 AT, 250 V AC, fusion lente, 5×20 mm), provenant du kit de fusibles

Outils nécessaires

Un tournevis à fente d'une taille comprise entre 3,3 mm et 5,5 mm convient

Préparatifs



AVERTISSEMENT—Décharge électrique

Des tensions élevées sont présentes dans l'appareil et peuvent causer des chocs électriques ou endommager l'appareil.

- Mettez l'appareil hors tension au moyen de son interrupteur d'alimentation principal. Déconnectez le cordon d'alimentation de la source d'alimentation et de l'appareil.
- Utilisez uniquement les fusibles du type et de courant nominal tels que spécifiés par Thermo Fisher Scientific pour l'appareil.
- N'utilisez pas des fusibles réparés et ne court-circuitez pas les portefusibles.

Procédez comme suit

Le porte-fusible est placé à proximité de l'interrupteur d'alimentation principal.



Illustration 43: Porte-fusible

N°	Description
1	Interrupteur d'alimentation principal (marche/arrêt)
2	Porte-fusible
3	Connecteur d'entrée d'alimentation

- 1. Retirez le porte-fusible à l'aide du tournevis.
- 2. Remplacez les deux fusibles par de nouveaux fusibles du type et du courant nominal indiqué. Remplacez toujours *les deux* fusibles.
- 3. Remettez en place le porte-fusible.
- 4. Rebranchez le cordon d'alimentation à la source d'alimentation et à l'appareil.
- 5. Mettez l'appareil sous tension au moyen de l'interrupteur d'alimentation principal.

7.9 Remplacement de la porte de l'appareil

À quel moment exécuter cette opération ?

La porte du compartiment à colonnes est défectueuse

Éléments nécessaires

Porte pour compartiment à colonnes

Préparatifs

Ouvrez le capot de la chambre à colonnes. Reportez-vous à la section Accès aux composants internes () page 47).

- 1. Ouvrez la porte à un angle de 90° de façon à aligner les deux paumelles du boîtier sur les rainures de la porte.
- 2. Pour décrocher la porte, soulevez-la.
- 3. Inclinez légèrement la porte vers l'extérieur du boîtier, puis décrochez-la.



Illustration 44: Articulation de la porte

N°	Description
1	Paumelle du boîtier
2	Rainure de la porte
3	Broche

- 4. Pour installer la porte, alignez les rainures de la porte sur les paumelles du boîtier.
- 5. Insérez la paumelle et le gond inférieurs dans la rainure de la porte.

- 6. Soulevez légèrement la porte pour insérer la paumelle et le gond supérieurs.
- Poussez la porte vers le bas pour la verrouiller en place. Vous pouvez fermer la porte seulement si elle est correctement installée.

7.10 Mise à jour du microprogramme de l'appareil

La description de cette section de rapporte au système de gestion de données chromatographiques de Chromeleon 7.

À quel moment exécuter cette opération ?

La mise à jour du firmware de l'appareil peut se révéler nécessaire, par exemple, après publication d'une nouvelle version de ce firmware qui ajoute des fonctions ou corrige les problèmes d'une version précédente.

Matériel nécessaire

Version du firmware ou de Chromeleon selon les besoins

NOTE Toute nouvelle version du firmware est incluse dans la dernière version révisée de Chromeleon. Le nouveau firmware *n'est pas* transféré automatiquement à l'appareil lorsque vous installez la version révisée de Chromeleon.

Préparatifs

- 1. Lisez les notes de publication fournies avec le microprogramme et/ ou la version de Chromeleon.
- 2. Connectez l'appareil au logiciel Chromeleon.
- 3. Arrêtez toutes les opérations sur l'instrument associé à l'appareil.
- 4. Patientez jusqu'à ce que l'instrument soit inactif.

- 1. Démarrer le programme Instrument Configuration Manager.
- Exécutez une mise à jour du firmware depuis l'onglet General de la boîte de dialogue de configuration de l'appareil. Pour obtenir plus de détails, consultez l'*Aide de Chromeleon*.
 La mise à jour du micrologiciel peut nécessiter plusieurs minutes.

AVIS

Une mise à niveau vers une version antérieure ou une mise à jour incomplète du micrologiciel peut donner lieu à une perte de fonctionnalité ou à un dysfonctionnement de l'appareil.

- N'interrompez pas la communication entre le logiciel Chromeleon et l'appareil pendant la procédure.
- Au début de la procédure de mise à jour, un message affiche la version du firmware installée sur l'appareil et la version à transférer depuis le logiciel Chromeleon. Si la version du firmware installée sur l'appareil est plus récente que la version affichée sous Chromeleon, annulez le téléchargement.
- Surveillez l'Audit Trail du programme Instrument Configuration Manager pour vérifier si la mise à jour du microprogramme a réussi ou échoué.
- 4. En fonction de la situation :

Situation	Action
Mise à jour du micrologiciel réussie	Une nouvelle qualification de l'appareil peut être nécessaire. Consultez les notes de mise à jour.
Échec de la mise à jour du micrologiciel	Éteignez puis rallumez l'appareil. Procédez à nouveau à la mise à jour du micrologiciel.
La mise à jour du micrologiciel échoue à plusieurs reprises	Prenez contact avec l'assistance technique de Thermo Fisher Scientific.

7.11 Déménagement ou expédition de l'appareil

Si vous souhaitez déménager l'appareil à un autre endroit, ou si vous devez l'expédier, vous devez tout d'abord le préparer selon les besoins. Suivez les instructions de cette section.

Respectez les consignes de sécurité suivantes :

AVIS

Pour lever ou transporter l'appareil, saisissez-le par les côtés. Ne déplacez et ne soulevez jamais l'appareil par la porte avant. Cela risque d'endommager les portes ou l'appareil.

Procédez comme suit

- 1. Préparez l'appareil pour le déménagement. Reportez-vous à la section Préparation de l'appareil au déménagement (▶ page 129).
- 2. Cette étape dépend des critères ci-après :
 - Pour déménager l'appareil vers un nouvel endroit, suivez les instructions de la section Déménagement de l'appareil (> page 130).
 - Pour expédier l'appareil, suivez les instructions sous Expédition de l'appareil (> page 131).

7.11.1 Préparation de l'appareil au déménagement

Pour préparer l'appareil au déménagement, suivez la procédure ciaprès :

- Exécutez la procédure d'arrêt de longue durée de l'appareil. Reportez-vous à la section Arrêt de longue durée (▶ page 97).
- 2. Mettez l'appareil hors tension au moyen de son interrupteur d'alimentation principal, puis débranchez le cordon d'alimentation.
- 3. Débranchez tous les câbles et raccordements fluidiques vers les autres appareils.
- 4. Obturez les orifices des vannes au moyen de bouchons de raccordement.
- Retirez l'appareil de refroidissement post-colonne (s'il est installé). Pour en savoir plus, reportez-vous à la section Remplacement de l'appareil de refroidissement post-colonne (VH-C10 uniquement) (▶ page 118).

- 6. Retirez l'appareil de préchauffage (s'il est installé) :
 - Pour retirer l'appareil de préchauffage actif, reportez-vous à la section Remplacement de l'appareil de préchauffage actif (> page 120).
 - Pour retirer l'appareil de préchauffage passif, reportez-vous à la section Remplacement de l'appareil de préchauffage passif (> page 122).
- 7. Retirez la colonne :
 - a) Débranchez les capillaires de l'entrée et de la sortie de la colonne.
 - b) Enlevez le porte-colonne de la cloison centrale à l'intérieur de la chambre à colonnes.
 - c) Retirez le porte-colonne du raccord de la colonne.
- Retirez le compartiment à colonnes du système Vanquish : Saisissez le compartiment à colonnes par ses côtés, soulevez-le légèrement puis sortez-le des encoches des barres de montage.

7.11.2 Déménagement de l'appareil

Préparatifs

Préparez l'appareil pour le déménagement. Reportez-vous à la section Préparation de l'appareil au déménagement (▶ page 129).

- 1. Respectez les consignes de manutention et de levage en toute sécurité de l'appareil.
- 2. Déménagez l'appareil vers son nouvel emplacement.
- Installez et mettez en place l'appareil dans le système empilé. Suivez les instructions de montage du système empilé du Manuel d'utilisation du système Vanquish.
- 4. Configurez l'appareil :
 - a) Connectez l'appareil et configurez les raccordements fluidiques (reportez-vous à la section Installation (> page 41)).
 - b) Préparez l'appareil à l'utilisation (reportez-vous à la section Préparation de l'appareil à l'utilisation (▶ page 88)).
- 5. Avant de lancer une analyse, laissez l'appareil s'équilibrer et assurezvous qu'il est prêt à fonctionner.

7.11.3 Expédition de l'appareil

Préparatifs

Préparez l'appareil pour le déménagement. Reportez-vous à la section Préparation de l'appareil au déménagement (▶ page 129).



ATTENTION—Contamination éventuelle

Des substances dangereuses peuvent contaminer l'instrument durant l'utilisation et peuvent causer des blessures corporelles au personnel d'entretien.

- Décontaminez toutes les pièces de l'instrument que vous souhaitez renvoyer pour réparation.
- Remplissez et signez le formulaire de santé et de sécurité. Thermo Fisher refuse d'accepter des instruments pour la réparation dont le formulaire de santé et de sécurité est manquant, incomplet ou non signé.

Procédez comme suit

1. Suivez les instructions de déballage de ce manuel dans le sens inverse.

Utilisez exclusivement le matériel d'emballage et la boîte de transport d'origine. Si la boîte de transport d'origine n'est pas disponible, des boîtes et du matériel d'emballage adéquats peuvent être commandés auprès du service commercial de Thermo Fisher Scientific.

2. Si vous devez retourner l'appareil à Thermo Fisher Scientific pour une réparation en atelier, contactez le service technique Thermo Fisher Scientific local pour savoir comment procéder.

Redémarrage de l'appareil après expédition

Pour installer l'appareil après l'expédition, suivez les instructions de montage du système empilé du *Manuel d'utilisation du système Vanquish*.

8 Dépannage

Ce chapitre est un guide pour résoudre les problèmes qui pourraient survenir pendant l'utilisation de l'appareil.

8.1 Généralités sur le dépannage

Les fonctions suivantes vous aident à identifier et à éliminer la source des problèmes qui peuvent survenir pendant l'utilisation de l'appareil.

NOTE Pour obtenir des renseignements sur les problèmes susceptibles de survenir pendant l'utilisation d'un système Vanquish, reportez-vous au *Manuel d'utilisation du système Vanquish*.

Si vous ne parvenez pas à résoudre un problème en suivant les instructions données dans cette section ou si vous rencontrez des problèmes non répertoriés, prenez contact avec l'assistance technique de Thermo Fisher Scientific pour obtenir de l'aide. Reportez-vous aux coordonnées figurant au début de ce manuel.

Afin de faciliter l'identification de l'appareil, munissez-vous du numéro de série et du nom technique lorsque vous communiquez avec Thermo Fisher Scientific.

Voyants d'état

La barre à LED d'état, située sur la façade avant de l'appareil, et la LED **STATUS**, située sur le panneau de commande intérieur, permettent un repérage visuel rapide de l'état opérationnel de l'appareil. Si le microprogramme détecte un problème, les voyants d'état émettent une lumière rouge.

Alarmes

Les fuites constituent un risque potentiel pour la sécurité. Aussi, si un capteur détecte une fuite, un signal sonore se déclenche afin de vous alerter, en plus du message affiché sur l'Instrument Audit Trail, et les voyants d'état deviennent rouges. Suivez les instructions de ce manuel pour trouver et éliminer la source de la fuite.

Messages de l'Instrument Audit Trail

Si le firmware de l'appareil détecte un problème, alors ce dernier est signalé au système de gestion de données chromatographiques.

Le système de gestion de données consigne tous les événements de la journée en cours relatifs à l'utilisation de l'instrument dans une Instrument Audit Trail. L'Instrument Audit Trail est nommée avec la date du jour, au format yyyymmdd. Par exemple, l'Instrument Audit Trail du 15 mai 2019 est nommée 20190515.

Les Instrument Audit Trails se trouvent dans le ePanel Set (Audit ePanel). De plus, les Audit Trails de chaque instrument sont disponibles dans le dossier de l'instrument concerné, dans la vue Chromeleon 7 Console Data.

Les messages de l'Instrument Audit Trail sont précédés d'une icône. Cette icône indique la gravité du problème. Pour connaître les causes possibles et les actions correctives, reportez-vous à la section Messages (> page 136).

Erreur du firmware

Lorsqu'une erreur du firmware survient pendant le fonctionnement du module, un journal des exceptions est créé et répertorie les processus lors de cette erreur du firmware. Le firmware envoie le journal des exceptions à l'Instrument Audit Trail lorsque le module est connecté au système de gestion des données chromatographiques.

Dans ce cas, respectez les consignes suivantes :

- Avant d'effacer le journal, envoyez l'Instrument Audit Trail à l'assistance technique en tant que fichier **.cmbx**.
- Pour effacer le journal des exceptions et reprendre le fonctionnement du module, exécutez la commande ExceptionLogClear.

Pour obtenir plus d'informations, consultez l'Aide de Chromeleon.

8.2 Messages

Ce tableau répertorie les messages les plus courants concernant l'appareil, ainsi que les procédures de dépannage.

Chaque message consiste en un code numérique accompagné d'un texte. Le code numérique est l'identifiant unique du problème ; le texte peut varier selon les modèles. Veuillez noter les points suivants :

- Afin de faciliter la recherche d'un message, le tableau répertorie les messages en les triant par code.
- Si vous ne trouvez pas le code recherché, vérifiez le texte du message. Les deux messages "Unexpected module behavior" et "Module malfunction detected" peuvent être attribués à des codes différents. Reportez-vous au début du tableau pour plus d'informations.

NOTE Si vous ne parvenez pas à résoudre le problème en suivant les instructions de ce manuel, ou si vous rencontrez un message non répertorié dans le tableau, notez-le en relevant son code et le texte et contactez-nous. Pour plus d'informations, reportez-vous à la section *Nous contacter* figurant au début de ce manuel.

Message et code	Description et action corrective
Unexpected module behavior. Code xx	Comportement inattendu du module. xx = code à deux, trois ou quatre chiffres. Lorsque le message s'affiche, notez son code, puis mettez le module hors tension. Patientez 5 secondes, puis remettez le module sous tension.
Module malfunction detected. Code xx	Détection d'un dysfonctionnement du module. xx = code à deux, trois ou quatre chiffres Lorsque le message s'affiche, notez son code. Mettez le module hors tension, puis prenez contact avec l'assistance technique.
Code 26 X tag path short circuit detected.	 Court-circuit chemin d'identificateur X détecté. (où X = 1 [appareil de préchauffage droit] ou 2 [appareil de préchauffage gauche]) 1. Débranchez le câble de l'appareil de préchauffage. 2. Si le message s'affiche de nouveau, prenez contact avec l'assistance technique.
Code 36 Download failed.	Échec du téléchargement. Le téléchargement du microprogramme a échoué. Relancez le téléchargement.

Message et code	Description et action corrective
Code 37 Download firmware mismatch.	Le microprogramme téléchargé ne correspond pas. Le téléchargement du microprogramme a échoué. Vérifiez que le fichier de microprogramme sélectionné est correct. Relancez le téléchargement.
Code 52 Module software incomplete. Download firmware (again).	Logiciel de module incomplet. Téléchargez le microprogramme (de nouveau). Le microprogramme est incomplet, par exemple parce que la communication entre le système de gestion de données chromatographiques et le module a été interrompue pendant le téléchargement du microprogramme. Relancez le téléchargement.
Code 84 Could not move X to its	Impossible de mettre X en la position souhaitée.
destination position.	(où X = vanne inférieure ou supérieure)
Code 85 Could not move X to its destination position.	La vanne de commutation indiquée est peut-être défectueuse. Remplacez la vanne de commutation, reportez-vous à la section Remplacement de la vanne de commutation pour colonne () page 113).
	Si le message d'erreur s'affiche de nouveau, prenez contact avec l'assistance technique.
Code 86 Unexpected X behavior.	Comportement inattendu de X.
Code 87 Unexpected X behavior.	(où X = vanne inférieure ou supérieure)
	La vanne de commutation indiquée est peut-être défectueuse. Remplacez la vanne de commutation, reportez-vous à la section Remplacement de la vanne de commutation pour colonne (▶ page 113). Si le message d'erreur s'affiche de nouveau, prenez contact avec l'assistance technique.
Code 89 Liquid leak sensor missing.	Capteur de fuite manquant.
	Prenez contact avec l'assistance technique de Thermo Fisher Scientific pour obtenir de l'aide. Notez que le module peut être utilisé même lorsque le capteur de fuites est défectueux ou absent, si la détection des fuites est désactivée.
Code 90 Download firmware mismatch –	Mauvaise correspondance de microprogramme – version invalide.
invalid version.	Vous avez essayé de télécharger un microprogramme incompatible avec un numéro de version antérieur à celui actuellement installé dans le module. Une mise à niveau vers une version antérieure du microprogramme peut donner lieu à une perte de fonctionnalité ou à un dysfonctionnement du module. Si nécessaire, relancez le téléchargement avec une version de microprogramme ultérieure à celle actuellement installée dans le module.
Code 91 Data changed during read.	Modification des données pendant la lecture.
	 Cela peut être une erreur temporaire. Mettez le module hors tension. Patientez 5 secondes, puis remettez le module sous tension.
	 Le microprogramme est peut-être défectueux. Mettez à jour le microprogramme.

Message et code	Description et action corrective
Code 118 USB Buffer Overflow.	Dépassement de tampon USB.
	Il s'agit d'un problème logiciel. Le module produit des données plus rapidement que l'ordinateur sur lequel le système de gestion de données chromatographiques lit les données.
	 Dans le système de gestion de données chromatographiques, déconnectez et reconnectez le module.
	 Si cela ne résout pas le problème, mettez à jour le microprogramme ou la version du système de gestion de données chromatographiques.
	3. Si le problème persiste : Un logiciel tiers installé dans l'ordinateur, comme les scanners antivirus, ou une mauvaise performance de l'ordinateur, peut également être à l'origine du problème. Contactez le service informatique sur site.
Code 120 System interlink request timed	Délai de requête System Interlink dépassé.
out.	La communication avec le module a échoué. Le module n'a pas répondu dans les temps.
	Pour le module pour lequel le message apparaît :
	1. Mettez le module sous tension si cela n'est pas encore fait.
	 Vérifiez les branchements System Interlink du module. Vérifiez que les deux extrémités de tous les câbles System Interlink sont branchées.
	3. Si le message s'affiche de nouveau, remplacez les câbles System Interlink.
Code 126 Device lost system interlink	L'appareil a perdu la connexion à System Interlink.
connection.	Un câble System Interlink peut être débranché du module ou défectueux.
	 Vérifiez les branchements System Interlink du module. Vérifiez que les deux extrémités de tous les câbles System Interlink sont branchées.
	2. Si le message s'affiche de nouveau, remplacez les câbles System Interlink.
Code 128 System interlink data errors.	Erreurs de données de System Interlink.
	Un câble System Interlink peut être débranché du module ou défectueux.
	 Vérifiez les branchements System Interlink du module. Vérifiez que les deux extrémités de tous les câbles System Interlink sont branchées.
	 Si le message s'affiche de nouveau, remplacez les câbles System Interlink.
Code 129 Device X not found on system	Appareil X introuvable sur System Interlink.
interlink.	X = numéro de série
	Le module avec ce numéro de série peut être débranché ou hors tension, ou une mise à jour du microprogramme peut être en cours.
	1. Vérifiez les branchements System Interlink du module.
	 Si le message s'affiche de nouveau, vérifiez l'état de fonctionnement du module.

Message et code	Description et action corrective
Code 131 System interlink bus not operable.	 Bus de System Interlink non opérationnel. Un câble System Interlink peut être débranché du module ou défectueux. 1. Vérifiez les branchements System Interlink du module. Vérifiez que les deux extrémités de tous les câbles System Interlink sont branchées. 2. Si le message s'affiche de nouveau, remplacez les câbles System
Code 136 Lock request rejected – already locked by X.	Requête de verrouillage rejetée – déjà verrouillé par X. X = identificateur du support de verrou, avec adresse USB liée au
	système de gestion de données de chromatographie ou adresse System Interlink liée au contrôleur du système ou à un module Le module est déjà verrouillé par un autre logiciel (contrôleur du système ou système de gestion de données chromatographiques). Attendez que le module soit libéré de son état verrouillé.
Code 137 Lock by X expired.	Verrouillage par X expiré.
	X = identificateur du support de verrou, avec adresse USB liée au système de gestion de données de chromatographie ou adresse System Interlink liée au contrôleur du système ou à un module
	Informez Thermo Fisher Scientific de l'événement. Aucune action supplémentaire requise.
Code 145 Lock holder X lost.	Support de verrouillage X perdu.
	X = identificateur du support de verrou, avec adresse USB liée au système de gestion de données de chromatographie ou adresse System Interlink liée au contrôleur du système ou à un module
	Le verrou a été libéré automatiquement puisque son support est disparu du System Interlink.
	Verifiez les branchements System Interlink du module.
Code 152 Assignment or command rejected – device locked.	Attribution ou commande rejetée – appareil verrouillé. Le module n'est pas prêt à accepter des commandes ou des attributions de propriété, puisqu'il est verrouillé par un autre logiciel (contrôleur du système ou système de gestion de données chromatographiques).
	Attendez jusqu'à ce que le module est libéré par le logiciel actuel (contrôleur du système ou système de gestion de données chromatographiques).
Code 2033 Column-ID X operation failed.	Défaillance de l'ID de colonne X.
	(où X = connecteur d'ID de colonne A, B, C ou D)
	Vérifiez d'abord que l'identificateur de colonne est correctement installé. Ensuite, reconnectez l'identificateur de colonne. Si le message s'affiche de nouveau, le connecteur d'ID de colonne peut être défectueux. Remplacez le connecteur d'ID de colonne.
Code 2034 Column-ID X operation failed.	Défaillance de l'ID de colonne X.
	(où X = connecteur d'ID de colonne A, B, C ou D)
	D'abord, formatez le système de fichiers de l'identificateur de colonne sous Chromeleon. Si le message s'affiche de nouveau, le connecteur d'ID de colonne peut être défectueux. Remplacez le connecteur d'ID de colonne.

Message et code	Description et action corrective
Code 2035 Column-ID X operation failed.	Défaillance de l'ID de colonne X. (où X = connecteur d'ID de colonne A, B, C ou D) Vérifiez d'abord que l'identificateur de colonne est correctement installé. Ensuite, reconnectez l'identificateur de colonne. Si le message s'affiche de nouveau, le connecteur d'ID de colonne peut être défectueux. Remplacez le connecteur d'ID de colonne. Si le message d'erreur s'affiche de nouveau, prenez contact avec l'assistance technique.
Code 2048 X leak detected - eliminate within approx. xxx seconds.	 Fuite de X détectée – Éliminez la fuite dans les xxx secondes suivantes environ. (où X = liquide) Le capteur de fuites a détecté un excès de liquide dans la chambre à colonnes. Supprimez la cause du problème dans le délai indiqué dans le message. Pour en savoir plus, reportez-vous à la section Résolution de fuites de liquide (▶ page 141).
Code 2049 X leak detected.	 Fuite de X détectée. (où X = liquide) Le capteur de fuites a détecté un excès de liquide dans la chambre à colonnes. Supprimez la cause du problème. Pour en savoir plus, reportez-vous à la section Résolution de fuites de liquide (▶ page 141).
Code 2070 Pre-heater X operation failed. Code 2071 Pre-heater X operation failed. Code 2072 Pre-heater X operation failed.	 Défaillance de l'appareil de préchauffage X. (où X = appareil de préchauffage gauche ou droit) Procédez comme suit : Vérifiez que l'appareil de préchauffage actif est correctement installé. Reconnectez l'appareil de préchauffage actif si nécessaire. Si le message s'affiche de nouveau, l'appareil de préchauffage actif est éventuellement défectueux. Remplacez-le. Reportez-vous à la section Remplacement de l'appareil de préchauffage actif (▶ page 120).
Code 2088 Module malfunction detected (X).	Module défectueux détecté (X). (où X = compartiment supérieur, compartiment inférieur, appareil de refroidissement post-colonne ou alimentation principale) Lorsque le message s'affiche, notez son code, puis mettez le module hors tension. Déconnectez tous les appareils de préchauffage actifs et mettez le module sous tension à nouveau. Si aucune erreur n'est signalée par le compartiment à colonnes par la suite, les appareils de préchauffage déconnectés sont défectueux. Remplacez-les. Reportez-vous à la section Remplacement de l'appareil de préchauffage actif (> page 120). Si le problème persiste, prenez contact avec l'assistance technique.

8.3 Résolution de fuites de liquide

À quel moment exécuter cette opération ?

Le capteur de fuites est humide. Le capteur de fuites signale une fuite.

Pièces et éléments complémentaires nécessaires

- Pièce de rechange, selon la nature du besoin
- Chiffon ou serviette en papier

Préparatifs

Lorsque vous essayez d'arrêter une fuite, respectez les consignes de sécurité et les règles générales de maintenance et d'entretien exposées à la section Maintenance et entretien (▶ page 101).

- Localisez l'origine de la fuite. Les fuites surviennent généralement aux raccords ; aussi, inspectez visuellement tous les composants et raccords du trajet d'écoulement.
- 2. Si nécessaire, serrez ou remplacez le raccord ou composant concerné.
- Au moyen d'un chiffon ou d'une serviette en papier, absorbez tout le liquide accumulé dans le bac de fuite et sous le capteur de fuites. Prenez garde à ne pas tordre le capteur.
- 4. Laissez le capteur s'équilibrer à la température ambiante pendant plusieurs minutes.
- 5. Si plus aucune fuite n'est signalée, vous pouvez reprendre l'utilisation de l'instrument.

9 Caractéristiques techniques

Ce chapitre répertorie les caractéristiques techniques matérielles et les caractéristiques de performance, et comprend notamment des informations sur les matériaux utilisés dans le trajet d'écoulement de l'appareil.

9.1 Caractéristiques de performance

L'appareil affiche les performances suivantes :

Туре	Caractéristique technique
Capacité d'accueil	2 loges à colonne
Plage de pression	VC-C10 : 2 – 70 MPa (20 – 700 bar, 290 – 10100 psi) VH-C10 : 5 – 151 MPa (50 – 1517 bar, 700 – 22002 psi)
Plage de température (chambre à colonnes)	VC-C10 : 5 à 85°C, max. 18°C en dessous de la température ambiante VH-C10 : 5 à 120°C, max. 18°C en dessous de la température ambiante
Exactitude de température (chambre à colonnes)	± 0,5°C (jusqu'à 80°C)
Stabilité de température (chambre à colonnes)	± 0,05°C
Précision de température (chambre à colonnes)	± 0,1°C
Délai de chauffage (chambre à colonnes)	De 20 à 50°C en moins de 15 minutes, ± 1°C De 25 à 40°C en moins de 5 minutes (généralement), ± 1°C
Délai de refroidissement (chambre à colonnes)	De 50 à 20°C en moins de 15 minutes, ± 1°C
Vannes de commutation pour colonne	Vanne 2 positions 6 ports Vanne multi-position 6 positions 7 ports
Communication	1 port USB (USB 2.0, connecteur de type "B")
System Interlink	2 ports System Interlink (connecteurs RJ45-8)
Commande	Chromeleon 7 L'appareil peut être également piloté à partir d'autres systèmes de gestion de données. Pour en savoir plus, veuillez prendre contact avec le service commercial de Thermo Fisher Scientific. Panneau de commande doté de 4 boutons permettant d'exécuter certaines fonctions directement depuis l'appareil.
Matériaux du trajet d'écoulement	Vanne : céramique/titane ; capillaires : MP35N ou acier inoxydable (pour obtenir des détails, reportez-vous aux caractéristiques techniques des appareils de préchauffage et de refroidissement post-colonne) AVIS Pour obtenir des informations sur la résistance chimique des matériaux, consultez la littérature technique.
Informations sur les solvants et les additifs	Reportez-vous à la section Informations sur les solvants et les additifs () page 25).
Fonctions de sécurité	Capteur de fuites
Fonctions de bonnes pratiques de laboratoire (BPL)	Les fonctions Predictive Performance vous permettent de programmer les procédures de maintenance sur la base des conditions réelles de fonctionnement et d'utilisation de l'appareil.
	Tous les paramètres système sont consignés dans la Chromeleon Audit Trail.
Appareils de préchauffage actifs

Туре	Caractéristique technique	
Plage de température	40°C à la limite de température supérieure de la chambre à colonnes	
	5°C au-dessus de la température ambiante et/ou de la température de l'éluant à l'entrée de l'appareil de préchauffage actif	
Matériaux dans le trajet d'écoulement	MP35N	
Biocompatibilité	Oui	
Performances de chauffage	80°C à 2,5 mL/min	
	120°C à 1 mL/min	
Exactitude de la température	± 2°C (jusqu'à 80°C)	
Stabilité de la température	± 0,5°C	
Précision de température	±1°C	
Volume chauffé	0,9 μL	
Volume total	0,1 x 380 mm :	0,1 x 610 mm :
	3 μL	6 μL

Appareils de préchauffage passifs

Туре	0,18 x 530 mm ; acier inoxydable	0,1 x 530 mm, MP35N	0,1 x 680 mm, MP35N
Plage de température	Identique à la chambre à colonnes	Identique à la chambre à colonnes	Identique à la chambre à colonnes
Biocompatibilité	Non	Oui	Oui
Volume chauffé	3 μL	1 μL	1 μL
Volume total	15 μL	5 μL	6 μL

Appareils de refroidissement post-colonne (VH-C10 uniquement)

Туре	Caractéristique technique	
Plage de température	40°C à 80°C	
Matériaux dans le trajet d'écoulement	MP35N	
Biocompatibilité	Oui	
Performances de refroidissement	Jusqu'à 50°C à 2 mL/min et avec le compartiment à une température de 120°C	
Exactitude de la température	± 2°C (jusqu'à 80°C)	
Stabilité de la température	± 0,5°C	
Précision de température	± 1°C	
Volume refroidi	1 μL	
Volume total	0,1 x 240 mm : 3 μL	0,1 x 590 mm : 6 μL

9.2 Caractéristiques physiques

L'appareil affiche les caractéristiques physiques suivantes :

Туре	Spécification
Type d'utilisation	Utilisation en intérieur uniquement
Plage de température d'utilisation	5 °C à 35 °C
Plage de température d'entreposage	-20 °C à 45 °C
Plage d'humidité d'utilisation	20 % à 80 % d'humidité relative, sans condensation
Plage d'humidité d'entreposage	60 % d'humidité relative maximum, sans condensation
Altitude d'utilisation	2 000 m au-dessus du niveau de la mer maximum
Degré de pollution	2
Exigences relatives à l'alimentation électrique	100 – 240 V CA, ± 10 % ; 50/60 Hz; max. 310 W / 330 VA
Catégorie de surtension	П
Niveau de pression acoustique	< 50 dB(A), généralement < 45 dB(A)
Dimensions (hauteur × largeur × profondeur)	70 x 11,1 x 42 cm
Poids	13 kg sans vannes de commutation pour colonne ; 14 kg avec deux vannes de commutation pour colonne installées

10 Accessoires, consommables et pièces de rechange

Ce chapitre répertorie les accessoires par défaut expédiés avec l'appareil et les accessoires disponibles en option. Ce chapitre fournit également des renseignements sur le réapprovisionnement en consommables et en pièces de rechange.

10.1 Généralités

L'appareil doit être utilisé exclusivement avec les pièces de rechange, composants supplémentaires, options et périphériques spécifiquement autorisés et certifiés par Thermo Fisher Scientific.

Les accessoires, consommables et pièces de rechange sont toujours conformes à la norme technique la plus récente. Les références des produits sont donc sujettes à modifications. Sauf indication contraire, les pièces les plus récentes sont compatibles avec les pièces qu'elles remplacent.

10.2 Kit d'expédition

Le kit d'expédition comprend les éléments répertoriés dans le tableau. Le contenu de ce kit est sujet à modifications et peut différer des informations de ce manuel. Consultez la liste incluse dans le kit afin d'obtenir les informations les plus récentes, à la réception de l'appareil.

Kit d'expédition

Élément	Quantité dans l'expédition
Pied avant (comportant l'orifice d'évacuation)	1
Pied arrière	1
Vis	4
Tournevis type Torx T20	1
 Kit de porte-colonnes (6 porte-colonnes), comprenant : Porte-colonne gauche permettant de placer la colonne à gauche (2 pièces) Porte-colonne droit permettant de placer la colonne à droite (2 pièces) Porte-colonne double permettant de placer deux colonnes à la même hauteur (2 pièces) 	2
Câble USB, type A à B, USB 2.0 ultra-rapide, 1 m	1
Kit d'installation standard, comprenant : • Vis (4 pièces) • Écrous rectangulaires en T (4 pièces) • Entretoises (4 pièces) • Pièce transversale (2 pièces)	1
ID de colonne (jeu de 2 connecteurs d'ID de colonne)	1
Câble System Interlink (RJ45), 0,5 m	1

Pour obtenir les informations de réapprovisionnement, reportez-vous à la section Consommables et pièces de rechange () page 151).

10.3 Accessoires en option

Élément	N° de référence
Appareil de préchauffage actif, 0,1 x 380 mm, MP35N	6732.0110
Appareil de préchauffage actif, 0,1 x 610 mm, MP35N	6732.0150
Kit de conversion Plusieurs kits de conversion peuvent être nécessaires pour assembler la configuration étendue avec jusqu'à trois compartiments à colonnes. Seuls les techniciens d'entretien certifiés par	Contactez le service commercial
Thermo Fisher Scientific sont habilités à installer la configuration étendue comprenant jusqu'à trois compartiments à colonnes ou à installer le compartiment à colonnes sur le côté gauche du système.	
Appareil de préchauffage passif, 0,1 mm x 530 mm, MP35N (VH-C)	6732.0174
Appareil de préchauffage passif, 0,1 mm x 680 mm, MP35N (VH-C)	6732.0184
Appareil de préchauffage passif, 0,18mm x 530 mm, acier inoxydable (VC-C)	6732.0170
Appareil de refroidissement post-colonne, 1 μL, 0,1 x 240 mm (VH-C10 uniquement)	6732.0510
Appareil de refroidissement post-colonne, 1 μL, 0,1 x 590 mm (VH-C10 uniquement)	6732.0520
Vanne 2 pos. 6 ports VC-C10 Le compartiment à colonnes peut accueillir une ou deux vannes de commutation pour colonne.	6230.1520
Vanne 6 pos. 7 ports VC-C10 Le compartiment à colonnes peut accueillir une ou deux vannes de commutation pour colonne.	6230.1530
Vanne 2 pos. 6 ports 150 MPa Bio VH-C/VC-C Le compartiment à colonnes peut accueillir une ou deux vannes de commutation pour colonne.	6036.2520
Vanne 6 pos. 7 ports 150 MPa Bio VH-C/VC-C Le compartiment à colonnes peut accueillir une ou deux vannes de commutation pour colonne.	6036.2530
Vanne 6 pos. 7 ports 70 MPa NP VC-C	6230.2530

10.4 Consommables et pièces de rechange

Fusibles et câbles

Description	N° de référence
Kit de fusibles Le kit comprend les fusibles adaptés aux modules du système Vanquish. Pour le compartiment à colonnes, utilisez uniquement des fusibles à fusion lente 4 AT (5 x 20 mm), 250 V AC.	6036.0002
Câble USB, de type A à type B, USB 2.0 ultra-rapide Longueur de câble : 0,5 m	6720.8910A
Câble USB, de type A à type B, USB 2.0 à haut débit Longueur de câble : 1 m	6035.9035A

Divers

Description	N° de référence
Portes pour compartiments à colonnes	6732.3890
Connecteurs d'ID de colonne	6732.0610
Capot de vanne de commutation pour colonne	6732.1843
Capteur à immersion PT100 Pour qualification opérationnelle et des performances du compartiment à colonnes	6705.0060
Matériaux d'emballage	6732.7004
Kit de qualification des performances Pour qualification opérationnelle et des performances du compartiment à colonnes, comprenant un thermomètre P700 et un capteur à immersion PT100. Note : pour continuer à utiliser un ancien kit de qualification des performances réf. 5705.0050A, il faut commander l'attache de capteur de température.	6732.0010
Rotor 6 pos. 7 ports NP VC-C	6230.2537
Stator 6 pos. 7 ports NP VC-C	6230.2535
Attache de capteur de température Pour qualification opérationnelle et des performances du compartiment à colonnes	6732.0009
Filtre en ligne Viper Pour retenir les particules (d'une taille supérieure à 0,5 μm) de la phase mobile	6036.1045
Raccord-union Viper	6040.2304
Pour les appareils de préchauffage actifs, passifs et les appareils de r post-colonne, reportez-vous à la section Accessoires en option () pa	refroidissement ge 150).

Cordons d'alimentation

Description	N° de référence
Cordon d'alimentation, Australie	6000.1060
Cordon d'alimentation, Chine	6000.1080
Cordon d'alimentation, Danemark	6000.1070
Cordon d'alimentation, UE	6000.1000
Cordon d'alimentation, Inde, AS	6000.1090
Cordon d'alimentation, Italie	6000.1040
Cordon d'alimentation, Japon	6000.1050
Cordon d'alimentation, Royaume-Uni	6000.1020
Cordon d'alimentation, États-Unis	6000.1001
Cordon d'alimentation, Suisse	6000.1030

11 Annexe

Ce chapitre comporte des renseignements complémentaires sur la conformité.

11.1 Conformité aux normes et directives

11.1.1 Déclarations de conformité

Déclaration de conformité CE

L'appareil satisfait aux exigences requises pour le marquage CE et respecte les exigences en vigueur.

Déclaration de conformité EAC

L'appareil satisfait aux exigences requises pour le marquage EAC et respecte les exigences en vigueur.

Conformité aux directives RoHS

Ce produit est conforme aux exigences des directives RoHS (Restrictions of Hazardous Substances) :

 Directive RoHS européenne
 Directive portant sur l'utilisation limitée de certaines substances dangereuses dans les équipements électriques et électroniques

Le marquage CE apposé sur l'appareil atteste de la conformité du produit à la directive.

 Réglementation RoHS chinoise Mesures de contrôle de la pollution provenant de produits d'information électroniques

Le cas échéant, un des logos suivants peut être apposé sur l'appareil :

Logo	Description
	Le logo vert est apposé sur les dispositifs qui ne comportent pas les substances dangereuses répertoriées par la réglementation.
	Le logo orange, qui encercle une valeur à un ou deux chiffres, est apposé sur les dispositifs qui comportent des substances dangereuses répertoriées par la réglementation. Le nombre indique la durée d'utilisation sans risques pour l'environnement (environment-friendly use period, EFUP) du dispositif. Pendant cette période, le dispositif (si employé conformément à son utilisation prévue) ne présente pas de risques graves pour la santé humaine ou l'environnement. Pour plus d'informations, consultez le site http:// www.thermofisher.com/us/en/home/technical-resources/rohs-

Déclaration de conformité UKCA

L'appareil satisfait aux exigences requises pour le marquage UKCA et respecte les exigences en vigueur.

Conformité à la norme UL/CSA 61010-1

L'étiquette du laboratoire NRTL apposée sur l'appareil (par exemple, le marquage cTUVus ou CSA) indique que l'instrument satisfait aux exigences des normes applicables.

11.1.2 Conformité à la directive DEEE

Ce produit doit être conforme à la directive de l'Union européenne relative aux déchets d'équipements électriques et électroniques (DEEE). Cela est signalé par le symbole suivant :



Illustration 45: Symbole DEEE

Thermo Fisher Scientific a conclu des contrats avec une ou plusieurs entreprises de collecte et de recyclage des déchets dans chaque État membre de l'Union européenne (UE), et ces entreprises sont tenues de collecter ou de recycler ce produit. Pour obtenir plus d'informations, veuillez contacter Thermo Fisher Scientific.

11.1.3 Conformité au règlement de la FCC

Cet équipement a été testé et déclaré conforme aux limites pour un appareil numérique de classe A, conformément à la partie 15 du règlement de la FCC (Commission fédérale des communications) aux États-Unis.

Ces spécifications sont destinées à fournir une protection raisonnable contre les interférences nuisibles, lorsque l'équipement est utilisé dans un environnement commercial. Cet appareil génère, utilise et peut émettre de l'énergie par radiofréquence et, s'il n'est pas installé et employé conformément aux instructions, peut causer des interférences nocives avec les communications radio. L'utilisation de cet appareil dans une zone d'habitation est susceptible de causer des interférences nocives ; le cas échéant, l'utilisateur est contraint de corriger les interférences à ses frais.

11.1.4 Historique des versions du manuel

Révision	Produits couverts
4.0	VC-C10-A-03, VH-C10-A-03
3.0	VC-C10-A-03, VH-C10-A-0x
2.0a	VH-C10-A-02
2.0	VH-C10-A-02
1.0	VH-C10-A

Ces instructions ont été rédigées en anglais (instructions originales). Les autres versions linguistiques sont des traductions des instructions originales en anglais.

Index

Α

accessoires	147
en option	150
kit d'expédition	149
additifs	25
informations	25
alimentation électrique 4	45, 46
appareil	
arrêt	96
optimisation	95
préparation à l'utilisation	88
redémarrage après un arrêt de longue du	urée
	99
appareil de préchauffage actif	95
installation	73
remplacement	120
thermostatisation	92
appareil de préchauffage passif	
Installation	61
remplacement	122
appareil de refroidissement post-colonne	
Installation	59
remplacement	118
thermostatisation	91
arret	96
courte durée	96
longue durée	97
arrêt de longue durée	ספ דם
audit trail	/ ۲ ۱۵۸
duuit ti dii	134

B

Barre à LED	84,	86,	134
-------------	-----	-----	-----

С

capillaires	
installation	69
Viper	70
capot de la chambre à colonnes	
fermeture	49
ouverture	47
capteur de fuites	91, 141
caractéristiques techniques	143
performance	144

nhucique	140
physique	140
curiometeon	30 124
dualt train	134
configuration de rapparen	00
Example 2 Configuration du module	00
Smart Shutdown	111
Smart Standby	90
Smart Startup	90
classe de sécurité	09
	19
connecter	71
thermostatisation	/ 1 01
concontration on chloruro	
condensation	20
configuration	. 40, 55
disposition du système	51
	۲C
matériel	00
Conformité aux normes LIL/CSA	JI 155
conformité aux normes ou conformité avec la réglementation	155
connecteurs	20
consignes	
installation	42
maintenance	103
utilisation	105 83
consignes de sécurité	00
cas d'urgence	24
cordon d'alimentation	46
entretien	103
équipement de protection	20
généralités	19
installation	42
maintenance	103
gualification du personnel	20
risques généraux	22
sécurité électrique	21
utilisation	83
consommables	147
cordon d'alimentation	. 46, 59

D

déballage	38
décontamination	109
DEEE	155
démarrage	88
dépannage	133
généralités	134
messages	136
déplacement	44
Désactivation de l'alarme	85
détection de fuites	34
dispositif d'évacuation	55
disposition du système	51

Ε

entretien	101
équilibrage	88
équilibrage du système	88
Equilibration Time	90
erreur du firmware	135
évacuation	55
ExceptionLogClear (Chromeleon)	135
exigences	
cordon d'alimentation	46
exigences relatives au lieu d'installation	45
alimentation électrique	45, 46
condensation	46, 59
expédition	129

F

Fast Cool	. 85, 91
FCC	155
Fonctions Predictive Performance	111
fuite	141
fuites de liquide	141
fusibles	123
fusibles de l'alimentation principale	123

G

gants	. 21

ID de colonne	35, 66, 90
identification de colonne	35, 66
installation	41

consignes de sécurité	42
exigences relatives au lieu d'installation	45
système	43
Instrument Audit Trail	134
Interlink	57

K

kit d'expédition 149

L

LED D'ÉTAT	134
livraison	40
lunettes de sécurité	21

Μ

maintenance	101, 108
consignes de sécurité	103
décontamination	109
fonctions Predictive Performance	111
fusibles	123
intervalle	108
introduction	102
mise à jour du micrologiciel	127
nettoyage	109
optimisation	111
règles générales	106
Marquage CE	154
Marquage cTUVus	155
Marquage EAC	154
Marquage RoHS	154
marquage UKCA	155
matériel compris	40
messages	136
messages d'avertissement	16
Messages de l'Audit Trail	136
mise à jour du micrologiciel	127
mise en marche	79
mise sous/hors tension	87
mode air calme	31, 92
mode air forcé	32, 92
mode de thermostatisation	31, 92
air calme	31, 92
air forcé	32, 92

Ν

nettoyage 10)9
--------------	----

0

Ρ

panneau de commande	
pièces de rechange	147
plage pH	
plaque signalétique	17
porte	
remplacement	125
retrait	102, 125
présentation (fonctionnelle)	29
principe de fonctionnement	

R

raccordement	
cordon d'alimentation 59	9
raccordements fluidiques 68	8
Ready Temp Delta 91, 93	3
réapprovisionnement 147	7
redémarrage de l'appareil après un arrêt de	
longue durée 99	9

S

signaux de température	92
Smart Shutdown	96
Smart Standby	96
Smart Startup	89
solvant	
concentration en chlorure	26
informations	25
plage pH	26
symboles de sécurité 16,	17
System Interlink	57
Système de raccord Viper	70

Т

tampons	25
concentration	25
informations	25
Temps d'équilibrage	93
transport	129
tubes	69

U

Universal Serial Bus	57
USB	57
utilisation	81
consignes de sécurité	83
éléments de commande	84
interruption	96
mise sous/hors tension	87
utilisation prévue	18

V

vanne de commutation pour colonne	112
capot de vanne	116
fuite	112
installation	63
raccordement	76
remplacement	113
vanne inférieure	85
vanne supérieure	85
vêtements de protection	21
voyant d'état	
Barre à LED 84, 86,	134
LED d'état 84, 86,	134
vue interne	33

www.thermofisher.com

Thermo Fisher Scientific Inc. 168 Third Avenue Waltham Massachusetts 02451 USA

