

thermoscientific

Vanquish

Système Neo VN-A10, VN-C10, VN-P10, VN-S10

Manuel d'utilisation

4822.5001-FR Révision 2.0 • Septembre 2024





Copyright © 2024 Thermo Fisher Scientific Inc. Tous droits réservés.

Traduction du manuel d'utilisation original

Les descriptions de matériel contenues dans ce manuel correspondent aux appareils : VN-A10-A-02, VN-C10-A-01, VN-P10-A-01, VN-S10-A-01.

Marques déposées

Sauf indication contraire, toutes les marques de commerce citées sont la propriété de Thermo Fisher Scientific Inc. et de ses filiales.

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.

PharMed est une marque de commerce de Saint-Gobain Performance Plastics.

Torx est une marque de commerce de Acument Intellectual Properties, LLC.

VICI est une marque de commerce de Valco Instruments Co. Inc et de VICI AG International.

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.

Nous contacter

Vous pouvez trouver des informations sur tous les différents services d'assistance de Thermo Fisher Scientific en naviguant sur le site Web suivant :

https://www.thermofisher.com/us/en/home/technical-resources/ contact-us.html

Table des matières

1	Uti	lisation du présent manuel 19	5
	1.1	Présentation du présent manuel 10	6
	1.2	Conventions des messages de sécurité1	7
	1.3	Avis spécifiques	8
	1.4	Conventions typographiques	9
	1.5	Documentation de référence 20	0
2	Séc	urité 21	1
	2.1	Symboles de sécurité et messages d'avertissement	2
		2.1.1 Symboles de sécurité et messages d'avertissement de ce manuel 22	2
		2.1.2 Respect des consignes du présent manuel 22	2
		2.1.3 Symboles de sécurité sur le système 22	3
		2.1.4 Plaques signalétiques 23	3
	2.2	Utilisation prévue 2!	5
	2.3	Précautions de sécurité 20	6
		2.3.1 Informations générales relatives à la sécurité 20	6
		2.3.2 Qualification du personnel	7
		2.3.3 Équipements de protection individuelle 21	7
		2.3.4 Consignes de sécurité électrique 28	8
		2.3.5 Risques résiduels généraux 29	9
		2.3.6 En cas d'urgence	1
	2.4	Conformité aux normes et directives	2
3	Pré	sentation du système 33	3
	3.1	Caractéristiques du système 34	4
	3.2	Description du système	7
	3.3	Socle du système	8
		3.3.1 Pièces du socle du système	8
		3.3.2 Verrouillage du socle du système	9
	3.4	Contrôleur du système 40	0
	3.5	Pompe	1
		3.5.1 Composants internes de la pompe 4	1
		3.5.2 Principe de fonctionnement de la pompe 44	2

	3.6	Passeu	ır d'échantillon	44
		3.6.1	Composants internes du passeur d'échantillon	44
		3.6.2	Principe de fonctionnement du passeur d'échantillon	45
		3.6.3	Attributions des orifices de la vanne du passeur d'échantillon	48
		3.6.4	Compartiment à échantillons	49
	3.7	Compa	artiment à colonnes	53
		3.7.1	Composants internes du compartiment à colonnes	53
		3.7.2	Principe de fonctionnement du compartiment à colonnes	54
		3.7.3	Vue d'ensemble de la vanne de commutation pour colonne	54
	3.8	Bac à s	solvants	56
	3.9	Détect	ion des fuites	57
	3.10	Dispos	itif de rinçage de l'aiguille (vue d'ensemble)	58
	3.11	Systèn	ne de rinçage des joints (vue d'ensemble)	60
	3.12	Mise s	ous et hors tension	62
	3.13	Utilisa	tion du système	64
	3.14	Purge	(vue d'ensemble)	65
4	Dák			67
4	Dep	anage	2	07
	4.1	Embal	lage endommagé, constatation d'un défaut à la réception	68
	4.2	Matér	iel fourni	69
	4.3	Déball	age du système	70
	4.4	Déball	age du compartiment à colonnes	72
	4.5	Déverr	rouillage du socle du système	73
5	Inst	allatio	on	75
	51	Consig	nes de sécurité pour l'installation	76
	5.1	Evigon	ass relatives an light d'installation	70
	5.2	5 2 1		70 78
		5.2.2	Alimentation électrique	79
		5.2.3	Cordon d'alimentation	80
		5.2.4	Condensation	81
		5.2.5	Conditions d'utilisation	81
	5.3	Installa	ation du système	83
	5.4	Ouver	ture et fermeture des portes avant	85

	5.5	Installa	ation du compartiment à colonnes	. 86
		5.5.1	Installation du compartiment à colonnes dans le système empilé	86
		5.5.2	Installation de la vanne de commutation Vanquish	88
		5.5.3	Installation de la vanne de commutation VICI	90
		5.5.4	Installation de l'appareil de préchauffage passif	92
		5.5.5	Installation des colonnes dans le compartiment à colonnes	94
	5.6	Retrait	t des protections de transport du passeur d'échantillon	96
		5.6.1	Retrait du couvercle isolant	. 97
		5.6.2	Installation du couvercle isolant	. 98
	5.7	Installa	ation du dispositif d'affichage	. 99
	5.8	Conne	xion des câbles	106
		5.8.1	Vue d'ensemble du connecteur du module du système	106
		5.8.2	Vue d'ensemble du connecteur du contrôleur du système	107
		5.8.3	Connexion de câbles USB	108
		5.8.4	Connexion des câbles du dispositif d'affichage	111
		5.8.5	Connexion de câbles System Interlink	113
		5.8.6	Connexion d'un câble LAN	115
		5.8.7	Connexion de câbles d'alimentation	115
		5.8.8	Diriger le câble de signal dans les serre-câbles	117
	5.9	Conne	xion d'un spectromètre de masse	118
	5.10	Verrou	uillage du socle du système	119
	5.11	Mise e	n place des raccordements fluidiques	120
		5.11.1	Passage des capillaires et des tubes dans le système	120
		5.11.2	Raccordement des capillaires, des raccords et des tubes	121
		5.11.3	Raccordement des conduites d'évacuation	124
	5.12	Rempl	issage des réservoirs de solvant	129
	5.13	Rempl	issage des réservoirs de rinçage de l'aiguille	131
	5.14	Rempl	issage du réservoir de liquide de rinçage des joints	134
	5.15	Premiè	ère mise en marche du système	136
	5.16	Intégra	ation dans un appareil LAN	138
	5.17	Prépar	ation du système pour l'utilisation initiale	139
6	Util	isatio	n	141
	6.1	Consig	nes de sécurité pendant l'utilisation	142
	6.2	Inform	ations sur les solvants et les additifs	144
		6.2.1	Compatibilité générale	144
		6.2.2	Plages pH autorisées	144

	6.2.3	Concentrations autorisées	145
	6.2.4	Utilisation de solvants et d'additifs	145
6.3	Consig	nes générales pour l'utilisation du système	147
6.4	Mise e	n marche du système	148
6.5	Éléme	nts de commande	149
	6.5.1	Panneaux de commande	149
	6.5.2	Barres à LED	153
6.6	Vue d'	ensemble de l'interface utilisateur Vanquish	155
	6.6.1	Informations sur le contrôleur du système	155
	6.6.2	Icônes d'état de l'état général du système	155
	6.6.3	Icônes de la barre de navigation et autres icônes	156
	6.6.4	État opérationnel du système	156
	6.6.5	Volet Accueil	157
	6.6.6	Volet Script	157
	6.6.7	Volet Paramètres	157
6.7	Réglag	e du dispositif d'affichage	159
	6.7.1	Positions possibles du dispositif d'affichage	159
	6.7.2	Déplacement du dispositif d'affichage de la position frontale à la posit	ion
		latérale	160
6.8	Prépai	er le système à l'utilisation	161
	6.8.1	Rinçage du système	161
	6.8.2	Equilibrage du système	163
	6.8.3	Chargement du carrousel	165
	6.8.3 6.8.4	Chargement du carrousel Recommandations sur les bouchons de flacons et les couvercles des	165
6.0	6.8.3 6.8.4	Chargement du carrousel Recommandations sur les bouchons de flacons et les couvercles des plaques à puits	165 167
6.9	6.8.3 6.8.4 Optim	Chargement du carrousel Recommandations sur les bouchons de flacons et les couvercles des plaques à puits isation des performances	165 167 169
6.9	6.8.3 6.8.4 Optim 6.9.1	Chargement du carrousel Recommandations sur les bouchons de flacons et les couvercles des plaques à puits isation des performances Interdépendance entre la pression de la pompe et le débit	165 167 169 169
6.9	6.8.3 6.8.4 Optim 6.9.1 6.9.2	Chargement du carrousel Recommandations sur les bouchons de flacons et les couvercles des plaques à puits isation des performances Interdépendance entre la pression de la pompe et le débit Ajout de solvants personnalisés au tableau d'étalonnage	165 167 169 169 170
6.9	6.8.3 6.8.4 Optim 6.9.1 6.9.2 6.9.3 6.0.4	Chargement du carrousel Recommandations sur les bouchons de flacons et les couvercles des plaques à puits isation des performances Interdépendance entre la pression de la pompe et le débit Ajout de solvants personnalisés au tableau d'étalonnage Changement de type de solvant	165 167 169 169 170 172
6.9	6.8.3 6.8.4 Optim 6.9.1 6.9.2 6.9.3 6.9.4 6.9.5	Chargement du carrousel Recommandations sur les bouchons de flacons et les couvercles des plaques à puits isation des performances Interdépendance entre la pression de la pompe et le débit Ajout de solvants personnalisés au tableau d'étalonnage Changement de type de solvant Changement du liquide de rinçage des joints	165 167 169 169 170 172 172
6.9	6.8.3 6.8.4 Optim 6.9.1 6.9.2 6.9.3 6.9.4 6.9.5 6.9.6	Chargement du carrousel Recommandations sur les bouchons de flacons et les couvercles des plaques à puits isation des performances Interdépendance entre la pression de la pompe et le débit Ajout de solvants personnalisés au tableau d'étalonnage Changement de type de solvant Changement du liquide de rinçage des joints Changement du liquide de rinçage de l'aiguille Réglage des décalages du captour de débit de la pompe	165 167 169 170 172 172 172
6.9	6.8.3 6.8.4 Optim 6.9.1 6.9.2 6.9.3 6.9.4 6.9.5 6.9.6 6.9.7	Chargement du carrousel Recommandations sur les bouchons de flacons et les couvercles des plaques à puits isation des performances Interdépendance entre la pression de la pompe et le débit Ajout de solvants personnalisés au tableau d'étalonnage Changement de type de solvant Changement du liquide de rinçage des joints Changement du liquide de rinçage de l'aiguille Réglage des décalages du capteur de débit de la pompe Bincage de l'aiguille	165 167 169 170 172 172 172 173 174
6.9	6.8.3 6.8.4 Optim 6.9.1 6.9.2 6.9.3 6.9.4 6.9.5 6.9.6 6.9.7 6.9.8	Chargement du carrousel Recommandations sur les bouchons de flacons et les couvercles des plaques à puits isation des performances Interdépendance entre la pression de la pompe et le débit Ajout de solvants personnalisés au tableau d'étalonnage Changement de type de solvant Changement du liquide de rinçage des joints Changement du liquide de rinçage de l'aiguille Réglage des décalages du capteur de débit de la pompe Rinçage de l'aiguille Thermostatage du compartiment du passeur d'échantillon	165 169 169 170 172 172 172 173 174
6.9	6.8.3 6.8.4 Optim 6.9.1 6.9.2 6.9.3 6.9.4 6.9.5 6.9.6 6.9.7 6.9.8 6.9.9	Chargement du carrousel Recommandations sur les bouchons de flacons et les couvercles des plaques à puits isation des performances Interdépendance entre la pression de la pompe et le débit Ajout de solvants personnalisés au tableau d'étalonnage Changement de type de solvant Changement du liquide de rinçage des joints Changement du liquide de rinçage de l'aiguille Réglage des décalages du capteur de débit de la pompe Rinçage de l'aiguille Thermostatage du compartiment du passeur d'échantillon	165 169 169 170 172 172 172 173 174 175
6.9	6.8.3 6.8.4 Optim 6.9.1 6.9.2 6.9.3 6.9.4 6.9.5 6.9.6 6.9.7 6.9.8 6.9.9 Arrêt	Chargement du carrousel Recommandations sur les bouchons de flacons et les couvercles des plaques à puits isation des performances Interdépendance entre la pression de la pompe et le débit Ajout de solvants personnalisés au tableau d'étalonnage Changement de type de solvant Changement du liquide de rinçage des joints Changement du liquide de rinçage de l'aiguille Réglage des décalages du capteur de débit de la pompe Rinçage de l'aiguille Thermostatage du compartiment du passeur d'échantillon Thermostatage du compartiment à colonnes	165 169 169 170 172 172 172 173 174 175 175
6.9	6.8.3 6.8.4 Optim 6.9.1 6.9.2 6.9.3 6.9.4 6.9.5 6.9.6 6.9.7 6.9.8 6.9.9 Arrêt o	Chargement du carrousel Recommandations sur les bouchons de flacons et les couvercles des plaques à puits isation des performances Interdépendance entre la pression de la pompe et le débit Ajout de solvants personnalisés au tableau d'étalonnage Changement de type de solvant Changement du liquide de rinçage des joints Changement du liquide de rinçage de l'aiguille Réglage des décalages du capteur de débit de la pompe Rinçage de l'aiguille Thermostatage du compartiment du passeur d'échantillon Thermostatage du compartiment à colonnes du système	165 167 169 170 172 172 172 173 174 175 175 177
6.9	6.8.3 6.8.4 Optim 6.9.1 6.9.2 6.9.3 6.9.4 6.9.5 6.9.6 6.9.7 6.9.8 6.9.9 Arrêt o 6.10.1 6.10.2	Chargement du carrousel Recommandations sur les bouchons de flacons et les couvercles des plaques à puits isation des performances Interdépendance entre la pression de la pompe et le débit Ajout de solvants personnalisés au tableau d'étalonnage Changement de type de solvant Changement du liquide de rinçage des joints Changement du liquide de rinçage de l'aiguille Réglage des décalages du capteur de débit de la pompe Rinçage de l'aiguille Thermostatage du compartiment du passeur d'échantillon du système Interruption du fonctionnement pour une courte durée Interruption du fonctionnement pour une longue période	165 169 169 170 172 172 172 173 174 175 175 177

7	Ma	inten	ance et entretien	179
	7.1	Prése	ntation de la maintenance et de l'entretien	. 180
	7.2	Consi	gnes de sécurité pour la maintenance et l'entretien	. 181
	7.3	Règle	s générales de maintenance et d'entretien	. 184
	7.4	Maint	enance de routine et maintenance préventive	. 186
		7.4.1	Plan de maintenance	. 186
		7.4.2	Rinçage de la pompe pour éviter toute contamination	. 189
		7.4.3	Rinçage de la pompe en cas de contamination persistante uniquemer	nt
				. 190
		7.4.4	Nettoyage ou décontamination d'un module du système ou du systèr	me . 191
	7.5	Maint	enance de la tête de pompe	. 195
		7.5.1	Vue d'ensemble des éléments de la tête de pompe	. 196
		7.5.2	Remplacement de la tête de pompe	. 197
		7.5.3	Remplacement des pistons	. 201
		7.5.4	Nettoyage du piston	. 203
		7.5.5	Remplacement de la bague de support	. 204
		7.5.6	Remplacement du joint du piston	. 209
		7.5.7	Remplacement du joint de rinçage des joints	. 216
		7.5.8	Remplacement des joints de la tête de pompe	. 220
		7.5.9	Entretien des clapets de non-retour de la tête de pompe	. 226
	7.6	Maint	enance du débitmètre	. 232
		7.6.1	Remplacement du filtre du débitmètre	. 232
		7.6.2	Remplacement de la conduite d'évacuation du débitmètre	. 234
	7.7	Entre	tien de la conduite de solvant	. 237
		7.7.1	Vidange des conduites de solvant	. 237
		7.7.2	Remplacement des conduites de solvant entre le réservoir et la pomp	be
		773	Remplacement des conduites de solvant entre la vanne d'arrêt et la t	. 237 Ato
		7.7.5	de pompe	. 242
		7.7.4	Remplacement de filtre de conduite de solvant	. 243
	7.8	Entret	tien de la boucle d'échantillonnage	. 245
		7.8.1	Composants de la boucle d'échantillonnage	. 246
		7.8.2	Retrait de la boucle d'échantillonnage	. 246
		7.8.3	Installation de la boucle d'échantillonnage	. 247
		7.8.4	Finalisation de la procédure de remplacement de la boucle	
			d'échantillonnage	. 249

7.9 Entr	etien de l'unité d'aiguille 250
7.9.1	Composant de l'unité d'aiguille et du siège de l'aiguille 250
7.9.2	Verrouillage de l'unité d'aiguille 251
7.9.3	Déverrouillage de l'unité d'aiguille 252
7.9.4	Remplacement du siège de l'aiguille et de l'unité de l'aiguille 254
7.10 Entr	etien de la tête du dispositif de mesure 263
7.10	1 Composant de la tête du dispositif de mesure 263
7.10	2 Retrait et raccordement des capillaires des têtes des dispositifs de mesure
7.10	3 Remplacement de la tête du dispositif de mesure
7.11 Entr	etien de la pompe d'évacuation 268
7.11	1 Test des tubes de la pompe d'évacuation 268
7.11	2 Remplacement du tube de la pompe d'évacuation 270
7.12 Entr	etien du dispositif de rinçage de l'aiguille 273
7.12	1 Remplacement des conduites de rinçage entre le réservoir et l'orifice d'admission de la pompe de rinçage de l'aiguille
7.12	2 Remplacement de la conduite de rinçage entre l'orifice de refoulement de la pompe de rinçage de l'aiguille et l'orifice de rinçage
7.12	3 Remplacement des conduites de rinçage entre les orifices d'admission du réservoir et du dégazeur
7.12	4 Remplacement des conduites de rinçage entre l'orifice de refoulement du dégazeur vers la vanne de sélection
7.12	5 Remplacement du clapet de non-retour du rinçage de l'aiguille
7.13 Entr	etien du système de rinçage des joints 282
7.13	1 Test d'étanchéité du système de rinçage des joints 282
7.13	2 Remplacement de la conduite de rinçage du réservoir de liquide de rinçage des joints
7.13	3 Remplacement de la conduite d'admission de rinçage des joints 288
7.13	4 Remplacement de la conduite d'admission de rinçage des joints de la
7.13	5 Remplacement du tube péristaltique dans la pompe de rinçage des joints
7.13	 6 Remplacement des conduites de rinçage des joints entre la pompe de rinçage des joints et le détecteur de rinçage des joints
7.13	7 Remplacement du détecteur de rinçage des joints 298
7.14 Insp	ection des vannes du passeur d'échantillon à la recherche de fuites 301
7.15 Entr	etien du compartiment à colonnes 303
7.15	1 Remplacement de colonnes 303
7.15	2 Inspection des vannes de commutation pour colonne à la recherche de fuites

		7.15.3	Remplacement de la vanne de commutation Vanquish	305
		7.15.4	Remplacement de la vanne de commutation VICI	308
		7.15.5	Remplacement du joint du stator et/ou rotor de la vanne de commuta VICI	tion 311
		7.15.6	Remplacement de l'appareil de préchauffage passif	313
	7.16	Contrô	eleur du système et maintenance du dispositif d'affichage	314
		7.16.1	Remplacement du contrôleur du système	314
		7.16.2	Remplacement de la carte SD	321
		7.16.3	Remplacement du dispositif d'affichage	323
	7.17	Rempla	acement du bac à solvants	326
	7.18	Rempla	acement de l'insertion de bouteille	329
		7.18.1	Retrait de l'insertion de bouteille	329
		7.18.2	Installation de l'insertion de bouteille	330
	7.19	Rempla	acement des fusibles de l'alimentation principale et des portes	332
		7.19.1	Remplacement des fusibles de l'alimentation principale	332
		7.19.2	Remplacement des portes	333
	7.20	Rempla	acement du module à glissières	336
		7.20.1	Préparation du retrait de la pompe	336
		7.20.2	Retrait du module à glissières de la pompe	337
		7.20.3	Préparation du retrait du passeur d'échantillon	338
		7.20.4	Retrait du module à glissières du passeur d'échantillon	340
		7.20.5	Préparation du retrait du compartiment à colonnes	341
		7.20.6	Retrait du module à glissières du compartiment à colonnes	342
		7.20.7	Retour du module à glissières	344
		7.20.8	Installation du module à glissières	345
	7.21	Prépar	ation du transport du système	347
	7.22	Fixatio	n des poignées de manutention	350
8	Dép	anna	ge 3	353
	8.1	Généra	alités sur le dépannage	354
	8.2	Messa	ges	356
		8.2.1	Messages communs	356
		8.2.2	Messages concernant la pompe	362
		8.2.3	Messages concernant le passeur d'échantillon	373
		8.2.4	Messages concernant le compartiment à colonnes	378

	8.3	Dépan	nage du système	. 379
		8.3.1	Pics	. 381
		8.3.2	Pression	. 386
		8.3.3	Variation des temps de rétention	. 388
		8.3.4	Ligne de base	. 389
		8.3.5	Temps de chargement de la colonne de piégeage prolongé	. 390
		8.3.6	Régulation de la température	. 391
		8.3.7	Alimentation et communications (modules du système)	. 392
		8.3.8	Problèmes en lien avec le contrôleur du système, le dispositif d'affich et l'interface utilisateur	age . 393
	8.4	Rechei	rche de fuites de liquide dans le système	. 395
9	Cara	actéris	stiques techniques	397
	9.1	Caract	éristiques physiques	. 398
		9.1.1	Spécifications environnementales	. 398
		9.1.2	Exigences relatives à l'alimentation électrique	. 398
		9.1.3	Dimensions	. 399
		9.1.4	Niveau de pression acoustique	. 399
		9.1.5	Poids	. 400
	9.2	Caract	éristiques de performance	. 401
		9.2.1	Caractéristiques de performance du système	. 401
		9.2.2	Caractéristiques de performance de la pompe	. 402
		9.2.3	Caractéristiques de performance du passeur d'échantillon	. 403
		9.2.4	Caractéristiques de performance du contrôleur du système	. 404
		9.2.5	Caractéristiques de performance du compartiment à colonnes	. 405
10	Acc	essoir	es, consommables et pièces de rechange	407
	10.1	Généra	alités	. 408
	10.2	Kit d'e	xpédition	. 409
		10.2.1	Système	. 409
		10.2.2	Compartiment à colonnes	. 412
	10.3	Access	oires en option	. 413
	10.4	Conso	mmables et pièces de rechange	. 417
		10.4.1	Capillaires, accessoires des capillaires et évacuation	. 417
		10.4.2	Pièces de rechange et outils du système	. 418
		10.4.3	Pièce de rechange du contrôleur du système	. 419
		10.4.4	Solvants et systèmes de rinçage	. 419
		10.4.5	Pièce de rechange de la pompe	. 421
		10.4.6	Pièces de rechange du passeur d'échantillon	. 423

10.4.7	Pièce de rechange du compartiment à colonnes 424
10.4.8	Câbles d'interface et cordons d'alimentation 425
11 Annexe	
11.1 Confo	rmité aux normes et directives 428
11.1.1	Déclarations de conformité 428
11.1.2	Conformité à la directive DEEE 429
11.1.3	Conformité au règlement de la FCC 429
11.1.4	Historique des versions du manuel 430
11.2 E/S nu	mériques (affectation des broches) 431
Index	

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 fournit les instructions d'installation, de mise en place, de démarrage, d'arrêt, de fonctionnement, de maintenance et de dépannage de votre système Vanquish Neo.

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 au système ou la perte de données.

Notes sur les illustrations

Les descriptions de ce manuel correspondent à une configuration standard du système Vanquish Neo.

- Les illustrations de ce manuel ne sont fournies qu'à des fins de compréhension élémentaire. Elles peuvent varier du modèle du système ou du composant.
- Aucune réclamation ne peut découler des illustrations de ce manuel.

1.2 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.

1.3 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 du système 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 du système.

1.4 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.

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* dans ce manuel sont employés selon le point de vue d'un observateur qui dirigerait son regard vers la façade avant du système.

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).

1.5 Documentation de référence

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

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 matériel

Pour obtenir des informations sur les bonnes pratiques et la configuration des applications prises en charge par votre système Vanquish Neo, consultez le *Guide de l'utilisateur du système Vanquish Neo*.

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 du système.

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 le système.

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 du système, afin de vous familiariser avec le système et ce manuel. Ce manuel comporte des informations importantes relatives à la sécurité des utilisateurs, ainsi qu'à l'utilisation et à la maintenance du système.
- Gardez toujours le manuel à proximité du système 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 le système

Ce tableau répertorie les symboles de sécurité qui apparaissent sur le système 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 du système.

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
0	Indique un bouton d'alimentation électrique qui permet de mettre sous tension ou hors tension le système.
(USC)	Indique un bouton d'alimentation électrique qui permet de mettre sous tension ou hors tension le contrôleur du système.
\sim	Indique un courant alternatif.
	Indique que la surface s'échauffe lors de l'utilisation. Ne touchez pas ces surfaces lorsqu'elles sont chaudes.
	Indique un danger potentiel de pincement. Veillez à ne pas introduire vos mains afin d'éviter de les blesser.
	Indique une borne de mise à la terre
=	Peut être utilisé pour relier un appareil à la terre, par exemple une colonne.

2.1.4 Plaques signalétiques

Les plaques signalétiques sont placées sur chaque module du système, à proximité des connexions électriques. Elles indiquent le numéro de série, le numéro de référence, le nom du module, le numéro de révision, et la tension nominale de la ligne électrique et des fusibles.

Étiquette du numéro de série du système

De plus, une étiquette du numéro de série du système est située du côté intérieur (en bas à gauche) du tiroir du socle du système. Cette étiquette indique le numéro de série du système, le numéro de référence, le nom du système et le numéro de révision.

Étiquette de type

Une étiquette supplémentaire située sur le bac de fuite de chaque module du système indique le nom du module, le numéro de série, le numéro de référence, ainsi que le numéro de révision. Afin de faciliter l'identification de l'appareil, gardez les informations de l'étiquette à votre portée lorsque vous communiquez avec Thermo Fisher Scientific.

Contrôleur du système

La plaque signalétique se trouve sur le panneau court du contrôleur du système. Elle indique le numéro de série et la tension nominale.

Dispositif d'affichage

La plaque signalétique se trouve à l'arrière du dispositif d'affichage. Elle indique le nom du module, le numéro de série, le numéro de référence, le numéro de révision (le cas échéant) et la tension nominale.

2.2 Utilisation prévue

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

Le système doit être utilisé uniquement par une personne qualifiée et dans un environnement de laboratoire.

Le système Vanquish Neo et ses modules doivent être utilisés comme équipement de laboratoire général (General Laboratory Equipment, GLE).

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 Neo 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 rayonnage 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 du système.



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

- Utilisez le système uniquement dans le cadre de ses caractéristiques techniques.
- Employez exclusivement les pièces de rechange, composants supplémentaires, options et périphériques spécifiquement autorisés et certifiés pour le système par Thermo Fisher Scientific.
- Ne réalisez que les procédures décrites dans le présent manuel d'utilisation et dans les documents relatifs au système. Suivez toutes les instructions étape par étape et utilisez les outils recommandés pour la procédure.
- N'ouvrez pas le boîtier du système et des autres composants, à moins que cela ne soit pas 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 du système. 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 l (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 du système.

Installation

Seules des personnes compétentes sont autorisées à mettre en service le système 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 autorisé par Thermo Fisher Scientific installe et met en place le système, il est de la responsabilité de l'installateur d'assurer la sécurité du système.

Utilisation



Fonctionnement général

Le système est destiné à une utilisation dans un environnement de laboratoire par du personnel formé et qualifié.

Tous les utilisateurs doivent connaître les dangers liés au système 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.

2.3.5 Risques résiduels généraux

Lors de l'utilisation du système, veuillez prendre en considération les risques généraux suivants :



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 le système 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 le système en présence de gaz ou de fumées inflammables.



AVERTISSEMENT—Vapeurs inflammables et dangereuses

Des vapeurs inflammables ou dangereuses peuvent s'échapper des récipients d'échantillon mal fermés contenant des échantillons inflammables ou volatils, et s'accumuler à l'intérieur du passeur d'échantillons. Cette accumulation peut présenter des risques pour la santé et la sécurité, et fausser les résultats.

- Utilisez uniquement des flacons ou des plaques à puits rendus étanches aux gaz au moyen de bouchons, de tapis d'étanchéité ou de rubans d'étanchéité. Consultez la liste actuelle des systèmes de fermeture autorisés par Thermo Fisher Scientific.
- Avant toute utilisation, inspectez les flacons afin d'y détecter d'éventuels fissures ou défauts. N'utilisez pas de flacons fissurés ou présentant des défauts.



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 les modules du système de l'alimentation électrique.

2.4 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 la livraison du système, ses modules respectent toutes les normes de compatibilité électromagnétique (CEM) et de sécurité applicables, telles que décrites dans le présent manuel.

Toute modification que vous apportez aux modules du système peut potentiellement annuler la conformité avec une ou plusieurs de ces normes de sécurité et de compatibilité électromagnétique. Les modifications des modules du système 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.

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

3 Présentation du système

Ce chapitre vous présente le système ainsi que ses principaux composants.

3.1 Caractéristiques du système

Le système présente les caractéristiques suivantes :

Caractéristiques principales

- Système pré-installé offrant une grande flexibilité au flux de travail
- Pression de fonctionnement jusqu'à 150 MPa
- Débit entre 100 nL/min et 100 μL/min
- Fonctions intégrées de démarrage et de purge automatisés
- Système de contrôle continu en arrière-plan pour une durée d'indisponibilité minimale
- Fonctions de diagnostic pour un dépannage rapide et un temps de disponibilité accru
- Interface utilisateur tactile, offrant une surveillance et un contrôle du système, ainsi qu'une assistance de maintenance

Caractéristiques spécifiques à la pompe

- Pompe binaire chaque voie étant munie d'un double piston de série en parallèle
- Conception de piston flottant pour compenser les petites tolérances mécaniques comprises dans les spécifications et ainsi améliorer la robustesse de la pompe
- Un débitmètre soutenant des débits entre 100 nL/min et 100 μL/ min aucun changement à apporter à l'appareil, à l'exception des capillaires du système
- Système de rinçage des joints actif pour une durée de vie prolongée des pistons et joints de piston
- Vannes de purge automatisées pour démarrage et diagnostic faciles
- Fonctions intégrées de démarrage et de purge automatisés
- Vannes d'arrêt automatiques intégrées empêchant le déversement de solvant et garantissant une bonne aspiration des solvants
- Dégazeur intégré pour les liquides de rinçage de l'aiguille du passeur d'échantillon

Caractéristiques spécifiques au passeur d'échantillon

- Principe d'injection avec division de flux à faible transfert interéchantillons présentant deux liquides de rinçage de l'aiguille (faible et fort)
- Rinçage de l'intérieur de l'aiguille, du siège de l'aiguille et de la boucle d'échantillonnage avec des liquides de rinçage faible et fort
- Rinçage de la surface extérieure de l'aiguille avec des liquides de rinçage faible et fort pour un transfert interéchantillons LC-MS minimal
- Volumes d'injection jusqu'à 100 μL avec une seule passe
- Volumes d'injections supérieures à 100 μL grâce à la fonction de prélèvements multiples pour la préconcentration de grands volumes
- Fonctions intégrées de démarrage et de purge automatisés
- Détection du fond pour permettre l'aspiration de petites quantités avec une perte minimale d'échantillon
- Régulation de la température pour la thermostatisation du compartiment à échantillons
- Lecteur de code-barres placé à l'intérieur du passeur d'échantillon identifiant les portoirs d'échantillons et les plaques à puits à l'aide de codes-barres Vanquish dédiés
- Prise en charge des portoirs d'échantillons et plaques à puits au format spécifié par la Society for Biomolecular Sciences (empreinte SBS)
- Nettoyage des joints de la tête du dispositif de mesure prolongeant la durée de vie des joints
- Fonctionnalité de chargement intégrée

Caractéristiques spécifiques au compartiment à colonnes

- Chambre à colonnes (34 x 8 x 4 cm, L x H x D) offrant de l'espace aux colonnes, aux colonnes de piégeage et aux capillaires dans différentes configurations de flux de travail
- Espace pour jusqu'à deux valves de commutation à l'intérieur de la chambre à colonnes pour les applications avancées
- La chambre à colonnes et les composants à l'intérieur de la chambre peuvent être thermostatés sur une grande plage de températures, avec une température minimale de 5°C au-dessus de la température ambiante et une température maximale de 80°C sans vannes ou de 60°C avec les vannes Vanquish Neo à faible dispersion

- Conception optimisée permettant une circulation d'air et une distribution de chaleur uniformes
- Temps d'échauffement rapide (de 35°C à 65°C ±1°C en moins de 12 minutes)
- Régulation de la température avancée offrant une stabilité de température de ±0,05°C
3.2 Description du système

Le système Vanquish est conçu pour réaliser des analyses de chromatographie liquide (LC) à faible débit et pour se connecter à un spectromètre de masse.

Le système Vanquish peut comprendre les modules suivants :



Illustration 1: Système empilé

N°	Description
1	Socle du système
2	Contrôleur du système
3	Pompe
4	Passeur d'échantillon
5	Compartiment à colonnes (en option)
6	Bac à solvants
7	Dispositif d'affichage

3.3 Socle du système

Le socle du système soutient la pompe, le passeur d'échantillon, le contrôleur et le compartiment à colonnes en option.

3.3.1 Pièces du socle du système

Le socle du système est doté des éléments suivants :

- des boutons d'alimentation pour mettre sous tension ou hors tension tous les modules et le contrôleur du système ;
- Un tiroir pour ranger les outils et les petites pièces du système ;
- Un orifice d'évacuation pour brancher une conduite d'évacuation du système ;
- Des verrous pour bloquer le système en place ou permettre de le déplacer.



Illustration 2: Socle du système (vues avant et arrière)

N°	Description
1	Bouton d'alimentation de Vanquish System Controller
	Permet la mise sous tension et hors tension du controleur du système.
2	Bouton d'alimentation du système Permet la mise sous tension / hors tension de tous les modules du système connectés au socle du système par le port System Interlink.
3	Tiroir à outils Permet le rangement de tous les outils nécessaires à l'installation ou à la maintenance du système Vanquish. Il suffit de le pousser pour l'ouvrir ou le fermer.
4	Réservé pour utilisation ultérieure
5	Orifice d'évacuation du système Permet le raccordement de la conduite d'évacuation du système.
6	Verrous, à gauche et à droite du socle du système Pour bloquer le système en place ou permettre de le déplacer.

N°	Description
7	Port System Interlink Permet le branchement des autres modules pour commander leur mise sous tension / hors tension à partir du socle du système Vanquish.
8	Port VSC Pour connecter le contrôleur système.

3.3.2 Verrouillage du socle du système

Le socle du système est doté d'un mécanisme de verrouillage. Avant de déplacer le système, par exemple, en vue de l'installation ou de la maintenance, vous devez déverrouiller le socle du système. Pour l'empêcher de se déplacer pendant l'utilisation, vous devez le verrouiller.

Vous pouvez déverrouiller et verrouiller le socle du système à l'aide des outils de verrouillage fournis.



Illustration 3: Outil de verrouillage du socle du système

3.4 Contrôleur du système

Les pièces du contrôleur du système sont illustrées ci-dessous.



Illustration 4: Socle du système avec bouton VSC et contrôleur du système

N°	Description
1	Bouton d'alimentation VSC
2	Agrafes-ressort
3	Emplacement pour carte mémoire flash (SD)
4	Réservé pour une utilisation future

3.5 Pompe

3.5.1 Composants internes de la pompe

Les composants de la pompe accessibles par l'utilisateur se situent directement derrière les portes avant.



Illustration 5: Composants internes de la pompe

N°	Description
1	Panneau de commande à boutons, pour la commande de pompe
2	Tête de pompe A (tête de pompe gauche)
3	Attaches de tube pour diriger les conduites de solvant vers les têtes de pompe
4	Tête de pompe B (tête de pompe droite)
5	Vannes d'arrêt
6	Capteur de fuites
7	Étiquette signalétique, indiquant le nom du module, le numéro de série, le numéro de référence et le numéro de révision
8	Orifice d'évacuation
9	Détecteur de rinçage des joints (détecteur de gouttes)
10	Dégazeur pour les liquides de rinçage de l'aiguille du passeur d'échantillon
11	Pompe de rinçage des joints
12	Gaine de tubulure munie de guides de tubulure
13	Débitmètre, avec :
13a	Entrée du débitmètre B
13b	Orifice de refoulement de la pompe (orifice de refoulement du débitmètre)
13c	Conduite d'évacuation du débitmètre (Purge Out)
13d	Entrée du débitmètre A



3.5.2 Principe de fonctionnement de la pompe

Illustration 6: Principe de fonctionnement

N°		Description
1		Conduites de solvant : A et B
2		Pompe à débit contrôlé avec deux blocs de pompe identiques (voies de solvant), A et B
	2a	Blocs de pompe A et B, chaque bloc de pompe muni de :
	2b	Cylindre de travail
	2c	Capteur de pression de fonctionnement
	2d	Cylindre d'équilibrage
	2e	Capteur de pression principale La pression principale est la pression dans la tête de la pompe.
3		Débitmètre, le cas échéant pour chaque voie, avec :
	3a	Conduite d'admission du débitmètre (capillaire depuis la tête de la pompe)
	3b	Filtre du débitmètre
	3c	Vanne de purge automatique avec orifice d'évacuation
	3d	Capteur de débit
	3e	Capillaire de restriction
	3f	Capteur de pression du système
	3g	Orifice de sortie du débitmètre avec pièce fluidique en T (orifice de sortie de la pompe)
	3h	Sortie d'évacuation
4		Capillaire entre système

La pompe est une pompe binaire munie d'un débitmètre pour contrôler le débit :

- Chaque bloc de pompe est muni d'un cylindre de travail et d'un cylindre d'équilibrage. Les cylindres sont connectés en série. Le solvant circule successivement dans les cylindres.
- Chaque bloc de pompe fournit un débit constant.
 Le cylindre de travail distribue le liquide et le cylindre d'équilibrage est rempli simultanément. Le cylindre d'équilibrage sert de réservoir et de refoulement tandis que le cylindre de travail effectue la course d'aspiration. Le cylindre d'équilibrage sert de réservoir et de refoulement tandis que le cylindre de travail effectue la course d'aspiration.
- Les capteurs du débitmètre mesurent et contrôlent les débits partiels des blocs de la pompe.
- Les débits partiels sont combinés dans la pièce fluidique en T à la sortie du débitmètre. Le débit combiné fournit au débitmètre la vitesse de débit sélectionnée et la composition du solvant.

3.6 Passeur d'échantillon

3.6.1 Composants internes du passeur d'échantillon

Les composants accessibles par l'utilisateur se situent directement derrière les portes avant.



Illustration 7: Vue interne du passeur d'échantillon du VN-A10

N°	Description
1	Compartiment à échantillons avec carrousel
2	Pompe d'évacuation
3	Bac de fuite avec :
3a	Panneau de commande avec voyants d'état
3b	Trou de guidage du capillaire de la pompe
3c	Étiquette signalétique, indiquant le nom du module, le numéro de série, le numéro de référence et le numéro de révision
3d	Pied d'évacuation
3e	Orifices d'évacuation
3f	Capteur de fuites
4	Pompe de rinçage de l'aiguille
5	Unité d'injection (reportez-vous à la page suivante pour obtenir une vue détaillée)
6	Couvercle isolant
7	Orifice d'admission de rinçage des joints
8	Guides pour capillaires vers le compartiment à colonnes



Unité d'injection du passeur d'échantillon

Illustration 8: VN-A10 Vue détaillée de l'unité d'injection sans couvercle isolant

N°	Description
1	Vis de verrouillage d'expédition pour l'unité d'aiguille
2	Unité d'aiguille avec pousseur de flacon
3	Unité de boucle d'échantillonnage
4	Siège de l'aiguille
5	Vanne d'injection
6	Vanne de solvant
7	Clapet de non-retour
8	Vanne de sélection
9	Tête du dispositif de mesure
10	Capteur de pression

3.6.2 Principe de fonctionnement du passeur d'échantillon

Le principe de fonctionnement du passeur d'échantillon repose sur le mode d'injection avec division de flux. Avec ce mode d'injection, l'aiguille fait partie intégrante de la boucle d'échantillonnage. La configuration décrite ici concerne le mode d'injection directe.

Prélèvement d'échantillon, vannes dans la position bypass

- Au début du processus d'injection d'échantillon, l'aiguille se déplace du siège de l'aiguille vers l'échantillon prédéfini placé dans le carrousel. Ensuite, l'aiguille descend dans le récipient d'échantillonnage.
- Le dispositif de mesure utilise l'aiguille pour aspirer l'échantillon dans le récipient d'échantillonnage et l'intégrer dans la boucle d'échantillonnage. Ensuite, l'aiguille est retirée du récipient d'échantillonnage.
- Afin d'éliminer tout échantillon résiduel de l'aiguille entre les injections, vous pouvez rincer l'extérieur de l'aiguille. Dans ce cas, l'aiguille se déplace vers l'orifice de rinçage et descend dans celui-ci. La pompe de rinçage de l'aiguille alimente l'orifice en liquide de rinçage afin de rincer la surface externe de l'aiguille et éliminer les déchets.
- L'aiguille revient se positionner dans le siège d'aiguille ce qui assure l'étanchéité de la pointe de l'aiguille sous la pression du système.

Injection de l'échantillon, vannes dans la position d'injection (voir la figure)

 La vanne d'injection bascule sur la position d'injection pour ouvrir le trajet d'écoulement permettant au solvant en provenance du débit de la pompe de s'écouler à travers la boucle d'échantillonnage, le siège de l'aiguille et l'aiguille, et de transporter l'échantillon à travers la vanne d'injection sur la colonne.

Démarrage du gradient, vannes en position de bypass

 Quand l'échantillon a été poussé complètement hors du passeur d'échantillon, la vanne d'injection revient à la position de bypass. Le gradient peut commencer à fournir le plus petit volume de retard des gradients possible.
 REMARQUE : si applicable, le passeur d'échantillon peut être configuré pour garder la vanne d'injection dans la position d'injection pour débuter le gradient immédiatement. Ainsi, la boucle d'échantillonnage complète est impliquée et le volume de retard des gradients sera plus grand.

Processus de rinçage, vannes dans la position de rinçage

Quand la vanne d'injection est dans la position de bypass, la vanne du solvant se place en position de rinçage et l'aiguille se déplace du siège de l'aiguille à l'orifice de rinçage et y descend. Dans cette position, l'extérieur et l'intérieur de l'aiguille seront rincés avec du liquide de rinçage de l'aiguille fort provenant de l'orifice de rinçage (7) et du solvant de rinçage faible fourni par le dégazeur en ligne placé dans la pompe (13).



Illustration 9: Principe de fonctionnement du passeur d'échantillon VN, position d'injection

N°	Description
1	Carrousel
2	Boucle d'échantillonnage
3	Aiguille
4	Siège de l'aiguille
5	Orifice de rinçage
6	Pompe de rinçage de l'aiguille
7	Liquides de rinçage de l'aiguille (fort et faible)
8	Vanne d'injection
9	Vanne de solvant
10	Évacuation des déchets
11	Capteur de pression
12	Dispositif de mesure
13	Solvants de rinçage (faible et fort) provenant du dégazeur placé dans la pompe
14	Vanne de sélection
15	Pompe
16	Colonne

3.6.3 Attributions des orifices de la vanne du passeur d'échantillon

Dans la configuration d'injection directe, les orifices de la vanne d'injection sont attribués comme suit :

Illustration 10: Vanne d'injection avec capillaires connectés

Orifice	Composant raccordé
1	Connecteur Viper
2	Vanne de solvant
3	Siège de l'aiguille
4	Colonne (si installé : compartiment à colonnes)
5	Boucle d'échantillonnage
6	Dispositif de mesure
7	Pompe



Pour l'injection directe et les flux de travail Retenir-et-Éluer, les orifices de la vanne de solvant sont attribués comme suit :



Orifice	Composant raccordé
1	Connecteur Viper
2	Déchets
3	Vanne d'injection
4	Connecteur Viper
5	Clapet de non-retour
6	Capteur de pression
7	Déchets

3.6.4 Compartiment à échantillons

Le compartiment à échantillons du passeur d'échantillon abrite un carrousel équipé de quatre segments à code-couleurs ; chacun de ces segments comporte un porte-flacon à trois positions pour des flacons de 10 ml.

Le passeur d'échantillon prend en charge des portoirs d'échantillons et plaques à puits au format spécifié par la Society for Biomolecular Sciences (format SBS).

Le compartiment à échantillons propose les fonctionnalités suivantes :

- Thermostatisation du compartiment à échantillons, voir Thermostatisation du compartiment à échantillons (> page 50).
- Identification et vérification du type de portoir, détection des segments vides et gestion de l'inventaire, reportez-vous à la section Identification du type de portoir (> page 51).

3.6.4.1 Thermostatisation du compartiment à échantillons

La régulation de la température peut être activée afin de permettre la thermostatisation du compartiment à échantillons. La circulation d'air permet de refroidir ou de réchauffer le compartiment à échantillons pour atteindre la température désirée, ce qui assure un équilibrage précis de la température de l'échantillon.

Pour obtenir des performances de thermostatisation optimales, maintenez les portes avant du passeur d'échantillon fermées pendant l'analyse de l'échantillon. N'ouvrez les portes avant que si cela se révèle nécessaire, par exemple pour charger les récipients d'échantillonnage dans le passeur d'échantillonnage.

La plage de température pour la thermostatisation du compartiment à échantillons est comprise entre +4°C et +40°C. Les échantillons peuvent être refroidis au maximum jusqu'à 23 K en dessous de la température ambiante.

En fonction de l'humidité ambiante, un liquide de condensation peut apparaître pendant le refroidissement des échantillons. Une pompe d'évacuation (pompe péristaltique) incorporée dans le passeur d'échantillon évacue activement le liquide de condensation du compartiment à échantillons.

Le passeur d'échantillon est muni d'un couvercle isolant couvrant la boucle d'échantillonnage. Ce couvercle protège les composants internes du taux d'humidité élevé et de la forte condensation lors de la thermostatisation du compartiment à échantillons. Si le couvercle n'est pas installé correctement, la thermostatisation ne pourra être activée ou s'arrêtera automatiquement.

3.6.4.2 Carrousel

Le carrousel comporte quatre segments à code couleurs : rouge (R), vert (G), bleu (B) et jaune (Y).

Chaque segment du carrousel compte les zones suivantes :

- Zone de positionnement pouvant accueillir un portoir d'échantillons ou une plaque à puits
- Porte-flacon supplémentaire comportant trois positions pour des flacons d'un volume de 10 ml maximum, tels que les flacons de réactifs
 Le porte-flacon supplémentaire est placé à la gauche de la zone

réservée au portoir d'échantillons ou à la plaque à puits dans chaque segment.



Illustration 12: Vue détaillée d'un segment du carrousel

N°	Description
1	Étiquette de code-barres « vide »
	Étiquette indiquant au lecteur de code-barres que le segment est vide
2	Code couleurs pour le segment (<i>ici</i> : R pour le segment rouge)
3	Zone de positionnement du segment
	Espace servant à positionner un portoir d'échantillons ou une plaque à puits
	Comporte des angles d'alignement et des points d'alignement pour aligner le portoir d'échantillons ou la plaque à puits
4	Porte-flacon supplémentaire
	Pour trois flacons d'un volume de 10 mL maximum
5	Code-couleurs pour le porte-flacon supplémentaire (<i>ici</i> : R1 par exemple pour le premier porte-flacon du segment rouge)

3.6.4.3 Identification du type de portoir

Un lecteur de codes-barres placé à l'intérieur du compartiment à échantillons permet d'identifier automatiquement le type de portoirs d'échantillons grâce aux plaques à puits sur lesquelles figure un codebarres 2D Vanquish indiquant le type de portoir.

Pendant le fonctionnement de l'appareil, le lecteur de codes-barres inventorie les portoirs présents en lisant le code-barres de type de portoir Vanquish, s'il est présent. Le lecteur de codes-barres identifie automatiquement le type de portoir et son orientation. Cette information est envoyée au système de gestion de données chromatographiques. Pour permettre l'identification du type de portoir d'échantillons ou de plaque à puits, utilisez des portoirs d'échantillons et plaques à puits munis de tels codes-barres.

Chaque segment contient une étiquette à code-barres qui informe le lecteur de codes-barres que le segment est vide et qu'aucun portoir d'échantillons ou plaque à puits n'est installé.



Illustration 13: Portoir d'échantillons avec code-barres pour l'identification du type de portoir

N°	Description
1	Position d'échantillon A1
2	Code-barres Vanquish pour le type de portoir

3.7 Compartiment à colonnes

3.7.1 Composants internes du compartiment à colonnes

Les composants du compartiment à colonnes accessibles par l'utilisateur se situent directement derrière les portes avant.



Illustration 14: Composants internes du compartiment à colonnes

N°	Description
1	Panneau de commande
2	Vanne de commutation pour colonne gauche (si installée ; ici : vanne Vanquish)
3	Orifices d'ID de colonnes
4	Ventilateur
5	Capteur de température
6	Guides pour capillaires
7	Vanne de commutation pour colonne droite (si installée ; ici : vanne Vanquish)
8	Étiquette signalétique, indiquant le nom du module, le numéro de série, le numéro de référence et le numéro de révision
9	Positions d'installation de l'appareil de préchauffage passif
10	Supports à colonnes

3.7.2 Principe de fonctionnement du compartiment à colonnes

Le critère fondamental d'un compartiment à colonnes destiné à des applications de chromatographie en phase liquide est la capacité à maintenir la température configurée avec le plus de précision possible. Grâce à la conception avancée du compartiment à colonnes, il est possible de stabiliser la température avec une précision de ±0,1°C.

Pour obtenir des résultats de mesure de bonne qualité, la stabilité de la température est plus importante que la précision de la température. Le compartiment à colonnes offre une stabilité de température de ±0,05°C.

Des éléments chauffants réchauffent la chambre à colonnes et les composants internes pour régler la température. La conception optimisée permet une circulation d'air uniforme.

Cette fonctionnalité garantit que :

- la température de la phase stationnaire reste constante sur toute la longueur de la colonne ;
- la colonne et l'éluant ont la même température pendant l'analyse ;
- la séparation analytique est effectuée à la température réglée.

3.7.3 Vue d'ensemble de la vanne de commutation pour colonne

Le compartiment à colonnes peut accueillir une ou deux vannes de commutation pour les applications avancées, y compris la commutation de colonnes. Vous pouvez installer les vannes de commutation dans les positions illustrées dans la section Composants internes du compartiment à colonnes () page 53).

Les vannes de commutation suivantes sont disponibles :

- vannes de commutation Vanquish ;
- vannes de commutation VICI[™]

Il est possible d'utiliser une configuration combinant ces deux types de vannes. Vous pouvez installer une vanne de commutation Vanquish partout où une vanne VICI n'est pas déjà installée.

Là où aucune vanne n'est installée, un couvercle ferme l'ouverture.

Vannes de commutation Vanquish

Les vannes de commutation Vanquish sont constituées d'une pièce. La commande de vanne est intégrée dans l'assemblage de vanne.



Illustration 15: Vanne de commutation Vanquish : vue de devant (gauche), vue de derrière (droite)

Pour installer une vanne Vanquish, reportez-vous à la section Installation de la vanne de commutation Vanquish (> page 88).

Pour remplacer une vanne Vanquish, reportez-vous à la section Remplacement de la vanne de commutation Vanquish (> page 305).

Vannes de commutation VICI

La vanne VICI est composée de deux parties, la capsule de vanne (n° 1) et la bague de verrouillage de vanne (n° 2). La bague de verrouillage de vanne connecte la capsule de vanne à la bride de la vanne (n° 3) de la commande de vanne. La commande de vanne n'est pas intégrée dans l'assemblage de vanne. Elle est situé dans le boîtier et le client n'y a pas accès.

Vous pouvez seulement installer ou remplacer la vanne VICI. Pour installer la commande de vanne VICI, prenez contact avec l'assistance technique locale de Thermo Fisher Scientific.





N°	Description
1	Capsule
2	Bague de verrouillage
3	Bride (à l'intérieur du boîtier)

Pour installer une vanne VICI, référez-vous à la section Installation de la vanne de commutation VICI () page 90).

Pour remplacer une vanne VICI, référez-vous à la section Remplacement de la vanne de commutation VICI (> page 308).

3.8 Bac à solvants

2a 2b 2c 2c

Le schéma montre les pièces du bac à solvants.

Illustration 17: Pièces du bac à solvants

N	°	Description
1		Bac à solvants
	1a	Devant aimanté Permet d'attacher un porte-étiquette magnétique sur le devant du bac à solvants, par exemple, pour indiquer un nom de système spécifique. Des étiquettes sont fournies dans le kit d'expédition du système.
2		Insertion de bouteille ¹ Dispose d'emplacements pour les réservoirs suivants :
	2a	5 emplacements pour les réservoirs de 1 L
	2b	2 emplacements pour les réservoirs de 2 L
	2c	4 emplacements pour les réservoirs de 0,25 L
1 b	¹ : l'insertion de bouteille peut être facilement retirée. Si vous retirez l'insertion de bouteille, utilisez de préférence uniquement les réservoirs de liquides livrés avec le	

bouteille, utilisez de préférence uniquement les réservoirs de liquides livrés avec le système. Si vous utilisez d'autres réservoirs de liquide, comme des réservoirs avec une plus grande capacité, assurez-vous que la pile du système est stable et ne s'incline pas. Si vous doutez de la sécurité, installez un kit de stabilisation de la pile (reportez-vous à la section Kit d'expédition () page 409)).

3.9 Détection des fuites

Comme les fuites constituent un problème potentiel de sécurité, les modules de pompe et de passeur d'échantillon sont munis de capteurs de fuites.

Ces capteurs de fuites détectent toute fuite dans le système provenant des raccordements fluidiques. Le liquide est recueilli dans les bacs de fuite, puis canalisé vers l'orifice d'évacuation qui permet d'évacuer le liquide par le dispositif d'évacuation.

Lorsque le capteur de fuites détecte une fuite, les voyants d'état du module en question deviennent rouges et un signal sonore se déclenche afin de vous alerter.

Si la fuite de la pompe n'est pas arrêtée dans les 3 minutes, la pompe interrompt le débit.

Consultez également

Recherche de fuites de liquide dans le système (> page 395)

3.10 Dispositif de rinçage de l'aiguille (vue d'ensemble)

Le dispositif de rinçage de l'aiguille du passeur d'échantillon permet de rincer l'intérieur et l'extérieur de l'aiguille afin d'en éliminer tout échantillon résiduel. Les cycles de rinçage peuvent être compris dans la séquence de prélèvement des échantillons ou réalisés manuellement.

Les composants de rinçage de l'aiguille sont pré-installés à l'intérieur du passeur d'échantillon lors de l'expédition.

Rinçage de l'extérieur de l'aiguille

Le dispositif de rinçage de l'extérieur de l'aiguille contient deux pompes de rinçage qui fournissent du liquide de rinçage fort ou faible.

L'extérieur de l'aiguille peut être rincé automatiquement après le prélèvement d'un échantillon ou manuellement.

Rinçage de l'intérieur de l'aiguille

Le rinçage de l'intérieur de l'aiguille comprend le rinçage de l'intérieur et de l'extérieur de l'aiguille. Pour garantir un faible transfert interéchantillons, le siège de l'aiguille et la boucle d'échantillonnage sont aussi rincés. En plus des deux pompes de rinçage de l'aiguille utilisées pour le rinçage extérieur, le rinçage intérieur utilise un liquide de rinçage de l'aiguille faible qui coule à travers un dégazeur en ligne situé dans la pompe.



Illustration 18: Composants de rinçage de l'aiguille

N°	Description
1	Conduite de rinçage étiquetée SWP (rouge) depuis le réservoir de liquide de rinçage fort vers l'orifice d'admission de la pompe de rinçage de l'aiguille, utilisée pour le rinçage fort de l'extérieur et de l'intérieur de l'aiguille.
	Conduite de rinçage étiquetée WWP (verte) depuis le réservoir de liquide de rinçage faible vers l'orifice d'admission de la pompe de rinçage de l'aiguille, utilisée pour le rinçage faible de l'extérieur de l'aiguille.
2	Conduite de rinçage étiquetée S (rouge) depuis le réservoir de liquide de rinçage fort vers l'orifice d'admission du dégazeur situé sur la pompe, utilisée pour les procédures de purge et d'arrêt de courte durée. Conduite de rinçage W (verte) depuis le réservoir de liquide de rinçage faible vers l'orifice d'admission du dégazeur situé sur la pompe, utilisée pour le rinçage faible de l'intérieur de l'aiguille et pour la purge.
3	Orifice de rinçage
4	Conduites de rinçage depuis l'orifice de refoulement de la pompe de rinçage de l'aiguille vers l'orifice de rinçage
5	Pompe de rinçage de l'aiguille
6	Conduites de rinçage depuis l'orifice de refoulement du dégazeur situé sur la pompe vers le passeur d'échantillon.
7	Conduite d'évacuation de rinçage de l'aiguille

3.11 Système de rinçage des joints (vue d'ensemble)

Le rinçage des joints de piston prolonge la durée de vie des pistons et des joints en empêchant la cristallisation et les dépôts de particules sur leurs surfaces et joints. Le trajet d'écoulement du système de rinçage des joints transite par la tête du dispositif de mesure située dans le passeur d'échantillon et par les têtes de pompe. Le système est constitué des pièces suivantes :

- Dans la pompe : pompe de rinçage des joints (pompe péristaltique), conduites de rinçage des joints et détecteur de rinçage des joints (détecteur de gouttes)
- Dans le passeur d'échantillon : conduites de rinçage des joints
- Réservoir de liquide de rinçage des joints

Lors de l'expédition, les composants de rinçage des joints et les conduites sont installés.



Illustration 19: Vue d'ensemble du système de rinçage des joints

N°	Description	
1	Orifice d'admission de rinçage des joints (orifice de droite)	
2	Réservoir de liquide de rinçage des joints entre le réservoir et l'orifice d'admission	
3	Conduite d'admission de liquide de rinçage entre l'orifice d'admission et la tête du dispositif de mesure	
4	Tête du dispositif de mesure	
5 + 6	Conduite d'admission de la pompe rinçage des joints	
7	Pompe de rinçage des joints (pompe péristaltique)	
8	 Conduites de rinçage des joints, dans la pompe (tube fin), avec : Conduite de rinçage des joints entre la pompe de rinçage des joints et la tête de pompe gauche Conduite de rinçage des joints sur la tête de pompe gauche Conduite de rinçage des joints entre les têtes de pompes Conduite de rinçage des joints sur la tête de pompe droite Conduite de rinçage des joints sur la tête de pompe droite Conduite d'évacuation de rinçage des joints entre la tête de pompe droite et le détecteur de rinçage des joints 	
9	Détecteur de rinçage des joints (détecteur de gouttes)	

Consultez également

Remplissage du réservoir de liquide de rinçage des joints
 () page 134)

3.12 Mise sous et hors tension

Boutons d'alimentation et interrupteurs d'alimentation principaux disponibles

L'image illustre les dispositifs d'alimentation du système.



Illustration 20: Boutons d'alimentation et interrupteurs d'alimentation principaux

N°	Description
1	Bouton d'alimentation du système
2	Bouton d'alimentation du contrôleur du système
3	Interrupteurs d'alimentation principaux

Bouton d'alimentation du système

Si les interrupteurs d'alimentation principaux des modules du système sont activés, vous pouvez utiliser le bouton d'alimentation du système pour mettre tous les modules du système sous ou hors tension simultanément (à l'exception du contrôleur du système) pour faciliter cette action. Respectez les consignes suivantes :

État	Description
Activé	 <i>Tous</i> les modules du système connectés au socle du système au moyen de câbles System Interlink sont mis sous tension en même temps. Il est impossible de mettre un module du système sous tension à l'aide du bouton d'alimentation du système si l'interrupteur d'alimentation principal du module du système est désactivé.
Désactivé	 <i>Tous</i> les modules du système connectés au socle du système au moyen de câbles System Interlink sont mis hors tension en même temps. Il ne suffit pas d'appuyer sur le bouton d'alimentation du système pour couper complètement l'alimentation électrique des modules du système. Pour mettre un module du système complètement hors tension, vous devez utiliser son interrupteur d'alimentation principal. Mettez un module hors tension au moyen de l'interrupteur d'alimentation principal lorsque vous en recevez l'instruction, par exemple avant des opérations de maintenance.

Bouton d'alimentation du contrôleur du système

Le bouton d'alimentation du contrôleur du système est un boutonpoussoir complètement indépendant du bouton d'alimentation du système. Si le contrôleur du système est mis sous tension, le dispositif d'affichage est automatiquement allumé. Le contrôleur du système redémarre automatiquement après une coupure de courant.

Pour mettre le contrôleur du système et le dispositif d'affichage (le cas échéant) hors tension, utilisez le bouton d'alimentation du contrôleur du système et attendez que la lumière LED (diode électroluminescente) d'état s'éteigne.

3.13 Utilisation du système

Maintenance et surveillance du système à l'aide de l'interface utilisateur Thermo Scientific Vanquish

Vous pouvez configurer le système et surveiller l'état du système à l'aide de l'interface utilisateur Thermo Scientific Vanquish. Cette interface vous permet de comprendre rapidement l'état du système, d'effectuer des procédures de maintenance et de résoudre des problèmes. Vous pouvez faire tout ceci avec ou sans connexion à un système de gestion de données chromatographiques.

Sauf mention contraire, l'ensemble des descriptions de logiciel présentes dans ce manuel se réfèrent à l'interface utilisateur Vanquish.

Pour une vue d'ensemble de l'interface utilisateur Vanquish, reportezvous à la section Vue d'ensemble de l'interface utilisateur Vanquish (> page 155).

Utilisation à partir d'un système de gestion de données chromatographiques

Pour obtenir des informations sur la connexion et le contrôle du système à l'aide du système de gestion de données chromatographiques, veuillez prendre contact avec le service technique Thermo Fisher Scientific local.

3.14 Purge (vue d'ensemble)

La purge consiste à rincer le système ou les composants du système pendant un certain temps à un débit élevé. Cette action est requise dans différentes situations et pour des raisons différentes. La situation et la raison déterminent le script à utiliser. Les tableaux contiennent l'information concernant les situations les plus courantes. Vous débutez les scripts à partir de la Vanquish User Interface. Suivez les instructions du script, le cas échéant.

Pendant l'installation, avant l'utilisation

Raison	Script	Paramètre supplémentaire
Lors de la préparation du système à l'utilisation initiale	A02 – Démarrer automatiquement	Aucun
Pour redémarrer le système après un arrêt de longue durée	A02 – Démarrer automatiquement	Aucun

Pendant l'utilisation

Raison	Script	Paramètre supplémentaire
Pour remplacer un réservoir de solvant ou un réservoir de liquide de rinçage de l'aiguille du passeur d'échantillon quand le niveau de liquide des réservoirs est trop bas pour la prochaine analyse (le même type de liquide)	B01 – Modifier les liquides / solvants	Sélectionner le solvant ou le liquide de rinçage de l'aiguille touché et Rafraîchir Uniquement = activé.
Une ou plusieurs conduites de solvant ou conduites de liquide de rinçage du passeur d'échantillon sont vides	B01 – Modifier les liquides / solvants	Sélectionner le solvant ou le liquide de rinçage de l'aiguille touché et Rafraîchir Uniquement = désactivé.
Pour changer pour un autre type de liquide (solvant ou liquide de rinçage de l'aiguille du passeur d'échantillon), par exemple, en changeant de méthode analytique	B01 – Modifier les liquides / solvants	Sélectionner le solvant ou le liquide de rinçage de l'aiguille touché et Rafraîchir Uniquement = désactivé.

Pendant la maintenance

Raison	Script	Paramètre supplémentaire
Maintenance de la pompe : Au début de la procédure de maintenance : pour éliminer les substances nocives À la fin de la procédure de maintenance : pour éliminer l'air qui pourrait être piégé dans le système en raison de la procédure	C02 – Purger la pompe	Éléments à purger = Ensemble de la pompe Intensité = Standard
Maintenance du passeur d'échantillon : Pour éliminer l'air qui pourrait être piégé dans le système en raison de la procédure	CO4 – Purger l'échantillonneur	Éléments à purger = Ensemble de l'échantillonneur Intensité = Intense

Pendant le dépannage

Raison	Script
Pour éliminer l'air qui pourrait être piégé dans le système et qui pourrait, par exemple, causer des problèmes de pulsation pendant l'utilisation, des niveaux de bruit élevés ou des analyses non reproductibles	Dépend de la cause potentielle. Reportez- vous à la section Dépannage du système (▶ page 379).

4 Déballage

Le présent chapitre décrit comment déballer le grand contenant d'expédition contenant le système empilé de base pré-installé. Pour de l'information sur comment déballer le compartiment à colonnes en option, reportez-vous à la section Déballage du compartiment à colonnes (> page 72). Le présent chapitre vous informe aussi sur les éléments compris dans la livraison.

4.1 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, après le déballage, inspectez le contenu afin d'y déceler d'éventuels signes d'une détérioration mécanique qui aurait pu survenir pendant le transport.

Si vous soupçonnez que le contenu 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.

4.2 Matériel fourni

Le matériel suivant est fourni :

- Système prémonté
- Dispositif d'affichage
- Manuel d'utilisation du système (téléchargeable à partir du site Web sur les manuels client)
- Cordons d'alimentation
- Kit d'expédition
 Pour obtenir des informations détaillées sur le contenu du kit, reportez-vous à Kit d'expédition (> page 409).

4.3 Déballage du système



ATTENTION—Charge lourde, système encombrant

Le système est trop lourd ou encombrant pour être soulevé en toute sécurité par une seule personne. Afin d'éviter toute blessure corporelle ou détérioration du système, veuillez respecter les consignes suivantes :

- Le maniement physique du système, notamment son soulèvement ou son déplacement, requiert la présence d'au moins deux personnes.
- Pour déplacer ou transporter le système, utilisez les poignées de manutention qui y sont fixées. Ne déplacez et ne soulevez jamais le système par les portes avant. Cela risque d'endommager les portes ou les modules du système. Vérifiez que les poignées sont bien fixées et ne peuvent pas glisser du système.

AVIS

Déplacer le système empilé Vanquish, même sur une courte distance sur la paillasse de laboratoire, peut endommager le socle du système et altérer son fonctionnement.

Avant de déplacer le système empilé, déverrouillez le socle du système à l'aide des outils de déverrouillage.

Outils nécessaires

Tournevis type Torx[™] T30

Procédez comme suit

- 1. Transportez le système jusqu'à l'emplacement où il doit être installé.
- 2. Retirez toutes les sangles de la boîte de transport.
- 3. Soulevez le couvercle en carton et retirez-le.
- 4. Retirez le kit d'expédition.
- 5. Soulevez la pièce en mousse supérieure et retirez-la.
- 6. Retirez le système de la boîte de transport :
 - a) Saisissez le système par les poignées de manutention.
 - b) Sortez le système de sa boîte de transport en le soulevant doucement et délicatement.
- 7. Placez le système sur une surface stable en laissant en place les films protecteurs recouvrant les surfaces du système.

- Pour déplacer légèrement le socle de système et mieux accéder aux pièces du système, déverrouillez le socle du système à l'aide des outils de déverrouillage du socle. Reportez-vous à la section Déverrouillage du socle du système (> page 73).
- 9. Retirez tout emballage supplémentaire.
- 10. Facultatif : Retirez les poignées de manutention :
 - a) Desserrez les vis de fixation de chaque poignée de manutention jusqu'à ce que les poignées de manutention puissent se déplacer dans le rail. Ne retirez pas complètement les vis des poignées de manutention.
 - b) Décrochez les poignées de manutention des rails, en les faisant glisser vers l'arrière du système.



Illustration 21: Poignées de manutention glissant du rail gauche

11. Fixez les vis des poignées de manutention.

NOTE Conservez la boîte de transport, les poignées de manutention accompagnées des vis de fixation, ainsi que tout le matériel d'emballage. Vous en aurez besoin si vous souhaitez transporter le système à un autre emplacement.

12. Retirez les films protecteurs de toutes les surfaces recouvertes.

4.4 Déballage du compartiment à colonnes

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, après le déballage, inspectez le contenu afin d'y déceler d'éventuels signes d'une détérioration mécanique qui aurait pu survenir pendant le transport.

Si vous soupçonnez que le contenu 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 poignées de manutention. Ne déplacez et ne soulevez jamais l'appareil par les portes avant. Cela risque d'endommager les portes ou l'appareil.

Procédez comme suit

- 1. Placez la boîte de transport par terre, puis ouvrez-la.
- 2. Retirez le kit d'expédition de la boîte de transport.
- 3. Retirez l'appareil de la boîte de transport ; pour ce faire, saisissez l'appareil par les poignées de manutention. Sortez l'appareil de sa boîte de transport en le soulevant doucement et délicatement.
- 4. Placez l'appareil sur une surface stable.
- Retirez tout emballage supplémentaire. Certaines surfaces, dont les portes de l'appareil, sont recouvertes d'un film protecteur lors du transport. Laissez tous les films protecteurs recouvrant l'appareil en place jusqu'à ce qu'il soit correctement positionné dans le système empilé.
- Au moyen des poignées de manutention, transportez l'appareil jusqu'au site d'installation s'il n'y est pas déjà, puis placez-le dans le système empilé. Reportez-vous à la section Installation du compartiment à colonnes dans le système empilé (> page 86).
- 7. Retirez les films protecteurs de toutes les surfaces recouvertes selon le cas.
4.5 Déverrouillage du socle du système

- 1. Introduisez un outil de verrouillage dans l'orifice de verrouillage d'un côté du socle du système.



- 2. Introduisez un outil de verrouillage dans l'orifice de verrouillage de l'autre côté du socle du système.
- Déverrouillez le socle du système en faisant tourner les outils de verrouillage comme illustré ci-dessous.
 Les outils de verrouillage ne peuvent pas être retirés du socle du système lorsqu'ils se trouvent en position « déverrouillé ».



Illustration 23: Déverrouillage du socle du système

- 4. Déplacez le système.
- Verrouillez le socle du système (reportez-vous à la section Verrouillage du socle du système (▶ page 119)).

5 Installation

Ce chapitre précise les exigences relatives au lieu d'installation et décrit comment mettre en place, installer et configurer le système.

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 26).



AVERTISSEMENT—Aiguille d'injection à pointe acérée

L'aiguille d'injection est dotée d'une pointe très acérée qui peut causer des blessures cutanées.

Pour éviter les blessures, ne touchez jamais la pointe de l'aiguille.



ATTENTION—Charge lourde, instrument encombrant

Les modules du système sont trop lourds ou encombrants pour être soulevés en toute sécurité par une seule personne. Afin d'éviter toute blessure corporelle ou détérioration des modules du système, veuillez respecter les consignes suivantes :

- Le maniement physique des modules, notamment son soulèvement ou son déplacement, requiert la présence d'au moins deux personnes.
- La présence de plusieurs personnes est requise en particulier pour placer les modules dans le système empilé ou pour les en retirer.
- Pour déplacer ou transporter les modules, utilisez les poignées de manutention qui les accompagnaient lors de la livraison. Ne déplacez et ne soulevez jamais les modules par les portes avant, car cela endommagerait les portes ou les modules. Vérifiez que les poignées sont correctement fixées et qu'elles ne peuvent pas glisser du module.



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

Déplacer le système empilé Vanquish, même sur une courte distance sur la paillasse de laboratoire, peut endommager le socle du système et altérer son fonctionnement.

Avant de déplacer le système empilé, déverrouillez le socle du système à l'aide des outils de déverrouillage.

Instructions pour l'extension de la pile système

Respectez les consignes suivantes :



AVERTISSEMENT—Risque de basculement du système empilé

Une configuration en une seule pile peut dépasser la hauteur maximale autorisée. Si la hauteur de la configuration du système est supérieure, le système empilé risque de basculer, ce qui peut provoquer des blessures.

Si la hauteur de votre configuration est supérieure, vous devez stabiliser la pile à l'aide d'une des mesures de sécurité suivantes :

- Utilisez un kit de stabilisation de la pile (reportez-vous au tableau ciaprès).
- Mettez le système en place sur deux piles distinctes.

Le tableau suivant indique quel kit de stabilisation de la pile est recommandé pour chaque hauteur de pile. Les hauteurs sont mesurées de la surface de la table jusqu'au rail du bac à solvants. En raison des tolérances de production et d'installation, la hauteur réelle de la pile peut être jusqu'à 2 cm supérieure à la limite. Cette tolérance est acceptable.

Hauteur de la pile	Kit de stabilisation de la pile requis
Jusqu'à 100 cm	Aucun Exemple : configuration avec une pompe, un passeur d'échantillon, un compartiment à colonnes et un détecteur à longueurs d'onde variables
Jusqu'à 123 cm	Un des kits suivants : • Kit de serrage de paillasse • Kit de stabilisation du système empilé • IonBench avec kit de montage empilé
Jusqu'à 139 cm	Kit de serrage de paillasse
139 cm et plus	Mettez le système en place sur deux piles distinctes

5.2 Exigences relatives au lieu d'installation

Un environnement d'utilisation approprié est indispensable pour assurer une performance optimale du système.

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

- Veillez à ce que le lieu d'installation dispose de suffisamment de bornes de connexion pour tous les appareils du système.
- Utilisez le système uniquement dans des conditions de laboratoire appropriées.
- Pour connaître les caractéristiques techniques, reportez-vous à Caractéristiques techniques (> page 397).

5.2.1 Paillasse

Le système Vanquish est conçu pour être posé sur une paillasse de laboratoire. Si la paillasse doit supporter un système Vanquish complet et éventuellement d'autres instruments, elle doit être capable de supporter le poids de tous les appareils et instruments, y compris les solvants.

Pour connaître les dimensions et les poids, reportez-vous à Caractéristiques techniques (> page 397).

Exigences relatives à la paillasse

Pièce	Exigences
Lieu d'installation de la paillasse	La paillasse doit être placée dans un endroit sûr et horizontal, sans vibrations.
Paillasse	 Robustesse Stable (non inclinable) Hauteur permettant un accès commode à l'intérieur de chaque appareil du système.
	 Le dessus de la paillasse doit être sec, propre et résistant aux produits chimiques.
	 Si la pile du système doit être stabilisée avec le kit de serrage de paillasse, le dessus de la paillasse doit remplir l'une des conditions suivantes :
	 Épaisseur : entre 15 et 50 mm ou
	 Matériau : suffisamment souple pour les vis à bois



Exigences relatives à l'encombrement du système

Illustration 24: Encombrement du système empilé (à gauche : vue de devant du système empilé, à droite : vue de dessus du système empilé)

N°	Description
1	Minimum 10 cm
2	Minimum 30 cm
3	Minimum 5 cm
4	Minimum 15 cm
5	Minimum 35 cm

5.2.2 Alimentation électrique

Les blocs d'alimentation des modules s'adaptent à un vaste éventail d'alimentations électriques et acceptent toutes les tensions secteur de la plage spécifiée pour le module.



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.

L'appareil peut subir de courtes coupures de tension si l'alimentation électrique de l'équipement est coupée pour le protéger d'un comportement indéterminé. Lorsque la tension d'alimentation revient, l'alimentation électrique redémarre automatiquement l'appareil et la connexion au système de données chromatographiques doit être rétablie.

5.2.3 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.

5.2.4 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.

5.2.5 Conditions d'utilisation

	Lors de l'utilisation du système, assurez-vous que le lieu d'installation respecte les exigences générales suivantes relatives aux conditions et à l'environnement d'utilisation. Pour connaître les exigences relatives à la température et à l'humidité ambiantes, reportez-vous à Caractéristiques techniques () page 397).
Température	
	Les fluctuations de température peuvent altérer les performances du système. Évitez les endroits présentant des variations importantes de température et soumis à de forts courants d'air. Par exemple, n'exposez pas directement le système à la lumière du soleil et ne le placez pas à proximité de sources de chaleur ou de froid, ou encore sous une conduite d'air.
Humidité	
	L'humidité relative de l'environnement de travail est essentielle pour un bon fonctionnement du système. Utilisez le système dans la plage d'humidité indiquée, et en l'absence de condensation.
	Si le taux d'humidité est trop élevé, de la condensation peut se former et endommager les composants électroniques du système. Si le taux d'humidité est trop faible, de l'électricité statique peut s'accumuler et se décharger, ce qui raccourcit la durée de vie des composants électroniques.

Ventilation

Assurez-vous que le lieu d'installation soit bien ventilé en permanence afin d'éviter les risques potentiels pour la santé et pour la sécurité pouvant résulter de la manipulation de substances liquides, volatiles ou gazeuses dangereuses.

Vibration et chocs mécaniques

Les vibrations et les chocs mécaniques peuvent altérer les performances du système. Par conséquent, le lieu d'installation ne doit pas être soumis à des vibrations. Évitez les chocs mécaniques et de placer le système à des endroits où d'autres instruments provoquent des vibrations.

Interférence électromagnétique

 Le système a été testé et déclaré conforme aux exigences d'immunité pour les environnements basiques et les environnements industriels électromagnétiques, conformément à la norme EN CEI 61326-1.

Les sources électromagnétiques de fortes interférences peuvent altérer les performances du système. Évitez l'emploi d'équipements générant de forts champs magnétiques ou électriques à proximité du système et utilisez uniquement des équipements de laboratoire certifiés. Ne branchez pas d'appareils pouvant perturber l'alimentation électrique sur le même réseau que le système.

 Le système a été testé et déclaré conforme aux limites d'émissions pour les équipements de classe A, conformément à la norme EN CEI 61326-1.



ATTENTION

Cet équipement n'est pas destiné à être utilisé au sein d'environnements résidentiels et il peut ne pas fournir une protection adéquate de la réception radio dans de tels environnements.

5.3 Installation du système

Un technicien d'entretien agréé par Thermo Fisher Scientific installe et configure le système Vanquish Neo, 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 Neo 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 le système, il convient de procéder comme suit.

Les étapes du diagramme sont les étapes d'installation de base. Suivez les instructions détaillées de chaque étape du diagramme dans la séquence indiquée.





Ouverture et fermeture des portes avant 5.4

Vous trouverez les composants accessibles par l'utilisateur et les raccordements fluidiques dans les appareils directement derrière les portes.

Ouverture des portes avant

- 1. Ouvrez la porte gauche.
- 2. Ouvrez la porte droite.

Mécanisme d'ouverture du La porte avant gauche du passeur d'échantillon est équipée d'un passeur d'échantillon mécanisme qui ouvre la porte automatiquement lorsque le passeur d'échantillon est hors tension afin de ventiler correctement le compartiment à échantillons. La porte s'ouvre aussi automatiquement en cas de coupure de courant ou si vous débranchez le cordon d'alimentation alors que le passeur d'échantillon est sous tension.





Fermeture des portes avant

- 1. Fermez la porte droite.
- 2. Fermez la porte gauche.

Mécanisme de fermeture du Lorsque le passeur d'échantillon est sous tension, la porte avant gauche passeur d'échantillon peut être complètement fermée.

5.5 Installation du compartiment à colonnes

5.5.1 Installation du compartiment à colonnes dans le système empilé

Le compartiment à colonnes est une pièce en option du système Vanquish. Les modules sont placés dans le système empilé, leur disposition dépendant de la configuration du système.

Outils nécessaires

- Tournevis type Torx T10
- Tournevis type Torx T20

Préparatifs

Lorsque le système est déjà utilisé :

- 1. Arrêtez le débit de la pompe.
- 2. Mettez l'alimentation électrique du système hors tension.
- 3. Déconnectez les conduites de liquide des réservoirs.
- 4. Retirez les réservoirs du bac à solvants.

Procédez comme suit

- 1. Desserrez les deux vis qui fixent le support en tôle sur le passeur d'échantillon à l'aide d'un tournevis (Torx T10).
- 2. Retirez le support en tôle.

NOTE Gardez le support et les vis. Vous pourriez en avoir besoin pour un transport ultérieur.



Illustration 26: Retrait du support en tôle (ici : passeur d'échantillon)

 Soulevez la partie avant du bac à solvants et retirez-le en tirant vers l'avant.

- 4. Soulevez le compartiment à colonnes par les poignées de manutention. Placez le compartiment à colonnes sur le passeur d'échantillon à environ 5 cm de l'extrémité des rails. Poussez le compartiment à colonnes vers l'arrière jusqu'à entendre le clic de mise en place.
- 5. Desserrez les vis qui maintiennent les poignées de manutention jusqu'à ce que vous puissiez les déplacer sur le rail, à l'aide du tournevis (Torx T20). Ne retirez pas complètement les vis des poignées de manutention.



Illustration 27: Poignées de manutention

N°	Composant
1	Poignées de manutention
2	Vis de fixation (une par poignée de manutention)

6. Décrochez les poignées de manutention des rails, en les faisant glisser vers l'arrière du compartiment à colonnes.



Illustration 28: Dépose de la poignée de manutention du rail gauche en la faisant glisser

- Placez le bac à solvants sur le compartiment à colonnes, à environ
 5 cm de l'extrémité des rails. Poussez le bac à solvants vers l'arrière jusqu'à ce qu'il s'enclenche en position.
- 8. Connectez les conduites de liquide aux réservoirs.
- 9. Connectez l'orifice d'évacuation sur le compartiment à colonnes au dispositif d'évacuation.
- 10. Connectez le cordon d'alimentation, la clé USB et les câbles du System Interlink au compartiment à colonnes.

Consultez également

- Connexion de câbles d'alimentation () page 115)
- Connexion de câbles USB (page 108)
- Connexion de câbles System Interlink () page 113)
- Passage des capillaires et des tubes dans le système () page 120)

5.5.2 Installation de la vanne de commutation Vanquish

Vous pouvez installer la vanne de commutation Vanquish à droite ou à gauche si une vanne VICI n'est pas déjà installée.

Éléments nécessaires

vanne de commutation Vanquish

Outils nécessaires

Tournevis type Torx T10

Préparatifs

- 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.

Procédez comme suit

- 1. Pour retirer le couvercle, vissez un bouchon de raccordement des orifices des vannes dans l'ouverture au centre du couvercle.
- 2. À l'aide du tournevis, desserrez les 4 vis sur le couvercle.
- 3. Retirez le couvercle en tirant sur le bouchon de raccordement.



Illustration 29: Retrait du couvercle



4. Retirez le ruban qui scelle le connecteur de câbles.

Illustration 30: Retrait du ruban

5. Retirez le câble du boîtier et connectez le câble au connecteur situé sur la face arrière de la vanne de commutation.



Illustration 31: Connexion du câble sur le connecteur

- 6. Enfoncez le câble dans le boîtier.
- 7. Enfoncez délicatement la vanne de commutation dans le boîtier.



Illustration 32: Installation de la vanne de commutation Vanquish

- 8. Vérifiez que la vanne de commutation est installée dans la bonne direction : l'orifice n° 3 doit être situé en haut.
- 9. Serrez les 4 vis de la vanne de commutation.



Illustration 33: 4 vis sur la vanne de commutation

- 10. Retirez les bouchons de raccordement qui couvrent les orifices des vannes.
- 11. Mettez le compartiment à colonnes sous tension.
- 12. Exécutez le script requis pour modifier votre flux de travail et suivez ses étapes pour raccorder les conduites de liquide.
- 13. Inspectez les vannes visuellement pour voir si elles fuient.
- Exécutez le script : D02 Tester le serrage du système (ce qui devrait être testé = le système). Ce script vérifie aussi si les vannes de commutation pour colonne présentent des fuites.

5.5.3 Installation de la vanne de commutation VICI

Vous ne pouvez installer la vanne VICI qu'aux emplacements où une commande de vanne VICI est installée. Pour installer la commande de vanne VICI, prenez contact avec l'assistance technique de Thermo Fisher Scientific locale. Pour une vue d'ensemble de la vanne de commutation VICI, reportez-vous à la section Vue d'ensemble de la vanne de commutation pour colonne () page 54).

Prérequis

La commande de la vanne VICI et le cadre d'adaptateur sont installés (effectué par l'ingénieur d'entretien).

Éléments nécessaires

• vanne de commutation VICI

Outils nécessaires

Tournevis type Torx T10

Procédez comme suit

- 1. Pour retirer le couvercle, vissez un bouchon de raccordement des orifices des vannes dans l'ouverture au centre du couvercle.
- 2. À l'aide du tournevis, desserrez les 4 vis sur le couvercle.
- 3. Retirez le couvercle en tirant sur le bouchon de raccordement.



Illustration 34: Retrait du couvercle

- 4. Ne retirez pas encore les bouchons de raccordement qui couvrent les orifices des vannes.
- 5. Alignez les deux broches situées à l'arrière de la bague de verrouillage de la vanne avec les ouvertures correspondantes sur la bride de la vanne.



Illustration 35: Alignement de la bague de verrouillage et de la bride

- 6. Pressez la capsule de la vanne sur la bride de la vanne.
- Pour aligner l'axe avec la capsule, appuyez sur le bouton VALVE LEFT ou le bouton VALVE RIGHT sur le panneau de commande et maintenez-le enfoncé, en fonction de la position de la vanne. La fonction de vanne de captage s'active et vous pouvez relâcher le bouton.
- 8. Quand l'axe est aligné, la capsule de vanne s'enclenche dans la bride de la vanne.

- 9. Retirez les bouchons de raccordement qui couvrent les orifices des vannes.
- 10. Connectez les conduites de liquide, selon les besoins de l'application.
- 11. Configurez la vanne de commutation VICI dans le système de gestion de données chromatographiques.
- 12. Inspectez les vannes visuellement pour voir si elles fuient.
- Exécutez le script : D02 Tester le serrage du système (ce qui devrait être testé = l'échantillonneur). Ce script vérifie aussi si les vannes de commutation pour colonne présentent des fuites.

5.5.4 Installation de l'appareil de préchauffage passif

L'appareil de préchauffage passif est une pièce en option. Vous pouvez l'utiliser pour porter l'éluent à la température de la colonne avant que l'éluent n'y pénètre.

Cet appareil s'utilise avec des débits de > 50 μ l/min, surtout avec des colonnes ID de 1 mm avec une gamme de débit entre 50 et 100 μ l/min.

Éléments nécessaires

- Appareil de préchauffage passif
- 2 vis

Outils nécessaires

Tournevis type Torx T10

Procédez comme suit

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

NOTE

L'appareil de préchauffage passif peut être installé dans n'importe lequel des deux 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.



Illustration 36: Positions d'installation de l'appareil de préchauffage passif

- 2. Alignez les trous de vis sur l'appareil de préchauffage passif avec ceux de l'emplacement de montage.
- 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.
- 4. Raccordez le capillaire de sortie de l'appareil de préchauffage passif à l'entrée de la colonne.
- Raccordez le capillaire d'entrée de l'appareil de préchauffage passif au passeur d'échantillon ou à la vanne de commutation pour colonne. En connectant l'appareil de préchauffage passif au passeur d'échantillon, dirigez le capillaire à travers le guide de capillaire le plus bas.



Illustration 37: Guide de capillaire pour diriger le capillaire au passeur d'échantillon

NOTE

La température de l'appareil de préchauffage passif n'est pas contrôlée. Il n'est donc pas nécessaire de configurer cet appareil dans le système de gestion de données chromatographiques.

5.5.5 Installation des colonnes dans le compartiment à colonnes

Huit supports à colonnes sont installés dans la chambre à colonnes. Attachez les colonnes sur les supports à l'aide des attaches pour colonnes du kit d'expédition. Vous pouvez installer les attaches pour colonnes sur les supports à colonnes dans n'importe quelle position.



Illustration 38: Positions possibles de l'attache pour colonne

Éléments nécessaires

- Colonne, selon les besoins de l'application
- Attaches pour colonne
- Coussins en mousse autoadhésifs

Procédez comme suit

1. Pour retirer la bague de la partie inférieure de l'attache pour colonne, appuyez légèrement sur la bague dans la direction des flèches.



Illustration 39: Attaches pour colonne

N°	Description
1	Bague
2	Coussin en mousse
3	Partie inférieure

- 2. Des coussins en mousse autoadhésifs sont expédiés avec les attaches pour colonne. Pour protéger la colonne, attachez un coussin à l'intérieur de chaque partie inférieure.
- 3. Attachez la partie inférieure de l'attache pour colonne au support de colonne dans la position voulue et tournez-la de 90° dans le sens des aiguilles d'une montre.
- 4. Raccordez les capillaires à l'entrée et à la sortie de la colonne.
- 5. Appuyez sur la colonne pour l'entrer dans la partie inférieure de l'attache pour colonne et réinstallez la bague.



Illustration 40: Réinstallation de la bague

6. Acheminez les capillaires par l'un des guides pour capillaires illustrés dans la figure.



Illustration 41: Guides pour capillaires

- Pour configurer les propriétés de la colonne, exécutez les scripts : A03 – Définir le type de colonne de séparation et A04 – Définir les caractéristiques de la colonne de séparation (et A05 – Définir les caractéristiques de la colonne de piégeage, le cas échéant).
- 8. Démarrez le débit de la pompe.

5.6 Retrait des protections de transport du passeur d'échantillon

Outils nécessaires

Tournevis plat

Procédez comme suit

 Retirez tous les verrous d'expédition ou capots de protection du passeur d'échantillon comme requis. Les verrous d'expédition à retirer sont marqués.



2. Ouvrez les portes du passeur d'échantillon.

Illustration 42: Ouverture des portes du passeur d'échantillon

Retirez le couvercle isolant (reportez-vous à la section Retrait du couvercle isolant (▶ page 97)).

Conservez-le, il devra être remis en place pour utiliser l'appareil.

AVIS

Si la vis de verrouillage d'expédition n'est pas correctement desserrée, le passeur d'échantillon peut être endommagé. Assurez-vous qu'elle est complètement desserrée avant de mettre en marche le passeur d'échantillon.

4. Desserrez la vis de verrouillage d'expédition qui se trouve au-dessus de l'unité d'aiguille. Avec vos doigts, tournez la vis imperdable dans le sens inverse des aiguilles d'une montre jusqu'à ce qu'elle reste suspendue librement dans le ressort. L'utilisation d'un tournevis plat peut faciliter la manœuvre.



Illustration 43: Desserrage de la vis de verrouillage d'expédition

- Installez le couvercle d'isolation (reportez-vous à la section Installation du couvercle isolant (▶ page 98)).
- 6. Retirez le rivet à pression du verrou d'expédition du carrousel et retirez le verrou.

5.6.1 Retrait du couvercle isolant

Procédez comme suit

- Saisissez le bas du couvercle et tirez légèrement dessus pour dégager la rainure dans le couvercle du support de fixation du couvercle isolant.
- 2. Tirez avec précaution sur le couvercle pour le désengager de la partie supérieure du cadre du boîtier.
- 3. Retirez le couvercle du support de fixation du couvercle isolant.



Illustration 44: Retrait du couvercle isolant

5.6.2 Installation du couvercle isolant

Éléments nécessaires

Couvercle isolant

Procédez comme suit

 Dirigez adéquatement le capillaire de la boucle d'échantillonnage à travers le guide désigné dans le support de montage du couvercle isolant.



Illustration 45: Couvercle isolant avec support de montage (à gauche) ; support de montage avec guide pour capillaires de la boucle d'échantillonnage (à droite)

- 2. Alignez le couvercle avec l'unité de l'aiguille et la boucle d'échantillonnage à l'intérieur du panneau avant.
- 3. Poussez soigneusement la rainure du couvercle sur le support de montage du couvercle isolant pour insérer le couvercle. Ne coincez pas le capillaire de la boucle d'échantillonnage avec le couvercle.

5.7 Installation du dispositif d'affichage

Vous pouvez fixer le dispositif d'affichage sur le côté gauche d'un module du système empilé (à condition que le module respecte les exigences figurant au paragraphe *Prérequis*). Pour faire pivoter le dispositif d'affichage, maintenez un dégagement de 10 cm vers la gauche.

Prérequis

Vous pouvez uniquement fixer le dispositif d'affichage sur les modules dont le boîtier comporte une encoche angulaire. Le dispositif d'affichage ne convient pas aux boîtiers comportant une encoche en U.



Illustration 46: Boîtiers avec les encoches (à gauche : angulaire, à droite : en U)

Préparatifs

Déballez le dispositif d'affichage.

NOTE Conservez la boîte de transport ainsi que l'ensemble du matériel d'emballage. Vous en aurez besoin si vous souhaitez transporter le dispositif d'affichage à un autre emplacement.

Éléments nécessaires

- Ensemble du dispositif d'affichage comprenant :
 - 3 vis Torx[™]
 - Bloc de fixation
 - Bloc coulissant



Illustration 47: Bloc coulissant et bloc de fixation

N°	Description
1	Bloc coulissant
2	Bloc de fixation

- Butée du dispositif d'affichage
- Cache du câble

Outils nécessaires

Tournevis type Torx T10

Procédez comme suit

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

1. Vérifiez si la position du câble du dispositif d'affichage correspond à l'illustration ci-dessous.





N°	Description
1	Bras pivotant
2	Câble du dispositif d'affichage

2. Si la position du câble du dispositif d'affichage ne correspond pas à l'illustration, faites tourner le dispositif d'affichage dans le sens horaire ou antihoraire au niveau du bras pivotant jusqu'à ce que le câble d'affichage soit guidé comme dans la figure ci-dessus.



Illustration 49: Rotation du dispositif d'affichage au niveau du bras pivotant

 Desserrez les 2 vis supérieures et dévissez la vis inférieure de l'ensemble du dispositif d'affichage afin de retirer le bloc de fixation.



Illustration 50: Vis concernées sur l'ensemble du dispositif d'affichage

4. En partant de l'arrière du rail du module, insérez le bloc coulissant du dispositif d'affichage dans le rail sur la gauche du module et tirez l'ensemble vers l'avant du système jusqu'à ce qu'il atteigne la butée.



Illustration 51: Insertion du dispositif d'affichage

N°	Description
1	Ensemble du dispositif d'affichage
2	Bloc coulissant
3	Rail

5. Fixez l'ensemble du dispositif d'affichage sans forcer à l'aide des 2 vis dans les deux trous du haut.



Illustration 52: Fixation du dispositif d'affichage

6. Localisez l'encoche à l'intérieur du panneau gauche du boîtier.



Illustration 53: Encoche

 Placez le bloc de fixation à hauteur de l'encoche, comme représenté. Vérifiez que le trou du bloc de fixation correspond à celui du boîtier.



Illustration 54: Positionnement du bloc de fixation

 Tout en maintenant le bloc de fixation avec la main droite, fixez-le sans forcer à l'aide d'une vis, main gauche à l'extérieur. Utilisez l'un des trous indiqués par un cercle dans l'image ci-dessous.



Illustration 55: Fixation du bloc de fixation

- 9. Vérifiez que le dispositif d'affichage est en position horizontale. Si ce n'est pas le cas, alignez-le en procédant comme suit :
 - a) Déplacez légèrement le dispositif d'affichage jusqu'à ce qu'il soit en position horizontale.
 - b) Serrez fermement les 3 vis.

10. Si le dispositif d'affichage est fixé au module supérieur, placez la butée du dispositif d'affichage au niveau du bac à solvants comme représenté sur la figure.



Illustration 56: Butée du dispositif d'affichage

5.8 Connexion des câbles

5.8.1 Vue d'ensemble du connecteur du module du système

Les modules sont équipés des connecteurs suivants pour le branchement de l'alimentation électrique et des câbles de signal :



Illustration 57: Connecteurs

N°	Description
1	Port System Interlink
	Permet de commander la mise sous tension / hors tension du système depuis le socle du système Vanquish et la communication de l'appareil
	Reportez-vous à la section Connexion de câbles System Interlink (> page 113).
2	Ports d'entrée et de sortie numériques (Dig I/O)
	Permettent l'échange de signaux numériques avec des instruments externes. Pour la connexion à un spectromètre de masse, utilisez le Dig I/O 2 sur le passeur d'échantillon.
	Reportez-vous à la section Connexion d'un spectromètre de masse (> page 118).
3	Concentrateur USB (Universal Serial Bus) (connecteur de type "A")
	Permet de connecter d'autres modules au système Vanquish
4	Port USB (connecteur de type "B")
	Permet de se connecter à d'autres modules du système Vanquish ou à l'ordinateur sur lequel est installé le système de gestion des données chromatographiques
	Reportez-vous à la section Connexion de câbles USB (> page 108).
5	Plaque signalétique (exemple) indiquant le numéro de série, le numéro de référence, le nom du module, le numéro de révision, et la tension nominale de la ligne électrique et des fusibles
6	Interrupteur d'alimentation principal (marche/arrêt)
7	Interrupteur d'alimentation principal
8	Connecteur d'entrée d'alimentation

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.

5.8.2 Vue d'ensemble du connecteur du contrôleur du système

Les connecteurs et emplacements suivants sont fournis sur le panneau court du contrôleur du système :



Illustration 58: Connecteur sur le panneau court

N°	Description
1	Plaque signalétique
	Indique le numéro de série et la tension nominale.
2	Connecteur Vanquish System Controller (VSC)
	Permet la connexion au socle du système.
3	Port System Interlink
	Permet le raccordement au socle du système et à un module dans le
	système HPLC
4	Port Universal Serial Bus (USB) (connecteur de type "A")
	Désactivé – Ne pas utiliser
5	Port Local Area Network (LAN)
	Permet la connexion au réseau du site.
6	Interrupteur d'alimentation principal (marche/arrêt)
7	Port d'entrée d'alimentation

Les connecteurs et emplacements suivants sont fournis sur le panneau long du contrôleur du système :





N°	Description
1	Port de sortie d'alimentation du dispositif d'affichage
2	Port de données du dispositif d'affichage
3	Ports System Interlink Désactivés
4	Port USB (connecteur de type "A") Désactivé – Ne pas utiliser
5	Étiquette indiquant le numéro de référence, le nom du module et le numéro de révision.
6	Emplacement pour carte mémoire flash (SD)

5.8.3 Connexion de câbles USB

Toutes les connexions USB requièrent un câble USB standard de type A vers B.



Illustration 60: Câble USB

L'extrémité plate et rectangulaire est le connecteur "A". L'extrémité plus petite et hexagonale est le connecteur "B".

NOTE

- La longueur de câble USB standard est limitée à 5 mètres. Les dispositifs USB ne doivent pas être plus éloignés de l'ordinateur ou du prochain concentrateur USB que 5 mètres.
- Après avoir branché les câbles USB et avant de mettre l'appareil sous tension pour la première fois, vérifiez que le logiciel de chromatographie est bien installé sur l'ordinateur. Quand l'appareil est mis sous tension, les pilotes USB nécessaires sont recherchés automatiquement et le système d'exploitation Windows[™] peut alors détecter l'appareil.


Illustration 61: Connexions de câble USB (ici : avec compartiment à colonnes installé)

N°	Description
-	Connexion USB :
1	Entre la pompe et l'ordinateur
2	Entre le passeur d'échantillon et la pompe
3	Entre le compartiment à colonnes (si installé) et la pompe

Procédez comme suit

- 1. Branchez le connecteur "B" du câble USB au port **USB** sur le passeur d'échantillon, la pompe et le compartiment à colonnes (si installé).
- 2. Branchez le connecteur "A" du câble USB de la pompe sur le concentrateur **USB** (2.0 ou ultérieur) de l'ordinateur.
- 3. Branchez le connecteur "A" du câble USB de chacun des autres modules sur le concentrateur **USB** de la pompe.

Pour contrôler les modules utilisant un concentrateur **USB** situé sur un autre module, le module muni du concentrateur doit être sous tension.

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 le système.

5.8.4 Connexion des câbles du dispositif d'affichage

Outils nécessaires

Ciseaux pour découper le cache du câble

Procédez comme suit

- 1. Branchez le câble d'alimentation du dispositif d'affichage dans le port **Display DC OUT 24V** du contrôleur du système.
- Insérez le câble de données du dispositif d'affichage dans le port
 Display Data du contrôleur du système et serrez à la main.
- 3. Acheminez les câbles de données et d'alimentation du dispositif d'affichage dans le rail.



Illustration 62: Acheminement des câbles du dispositif d'affichage

4. Découpez le cache du câble à une longueur de 34 cm à l'aide des ciseaux.



5. Enfoncez fermement le cache du câble dans le rail.

Illustration 63: Montage du cache du câble

6. Fixez la partie verticale du trajet du câble à l'aide des serre-câbles du boîtier.



Illustration 64: Fixation des câbles à l'aide des serre-câbles

- 7. Replacez le système empilé dans la position d'utilisation.
- Verrouillez le socle du système à l'aide des outils de verrouillage (reportez-vous à la section Verrouillage du socle du système (> page 119)).

5.8.5 Connexion de câbles System Interlink

Les connexions System Interlink nécessitent l'utilisation d'un câble spécial (câble de catégorie 6).



Illustration 65: Câble System Interlink

Les connexions System Interlink entre les modules sont pré-installées, à l'exception du compartiment à colonnes en option.



Illustration 66: Connexions de câbles System Interlink

N°	Description
-	Connexion System Interlink :
1	Entre le socle du système et le contrôleur du système
2	Entre le contrôleur du système et la pompe
3	Entre la pompe et le passeur d'échantillon
4	Entre le passeur d'échantillon et le compartiment à colonnes (si installé)

Si un compartiment à colonnes est installé (en option)

- 1. Retirez le bouchon qui recouvre le port **System Interlink** de droite du passeur d'échantillon.
- Branchez l'une des extrémités du câble System Interlink sur le port System Interlink du compartiment à colonnes et l'autre sur le port System Interlink de droite du passeur d'échantillon (comme représenté dans la figure).

5.8.6 Connexion d'un câble LAN

Éléments nécessaires

- Câble LAN
- Carte réseau

Procédez comme suit

AVIS—Endommagement du module Vanquish

Les ports **System Interlink** et **LAN** utilisent le même type de connecteurs. Le raccordement d'un câble LAN, par exemple, au port **System Interlink** d'un module Vanquish, peut endommager ce module et le contrôleur du système.

Branchez uniquement les câbles sur les ports correspondants.

- 1. Branchez un câble LAN sur le port LAN du contrôleur du système.
- 2. Branchez l'autre connecteur du câble LAN sur le port LAN de carte réseau de l'ordinateur sur lequel est installé le système de gestion de données chromatographiques.

5.8.7 Connexion de câbles d'alimentation

La présente section contient de l'information détaillée sur comment connecter les modules à la source d'alimentation. Ne mettez pas encore le système ni les modules sous tension.

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.

Procédez comme suit

- 1. Vérifiez que l'interrupteur d'alimentation de chacun des appareils est positionné sur OFF (hors tension).
- 2. Branchez le cordon d'alimentation au connecteur d'alimentation sur chaque appareil.
- 3. Branchez les extrémités libres des cordons d'alimentation à une source d'alimentation appropriée.
- Insérez les câbles de signal et d'alimentation dans les serre-câbles (reportez-vous à Diriger le câble de signal dans les serre-câbles (▶ page 117)).
- 5. Si nécessaire, déplacez le système vers son emplacement définitif.
- Verrouillez en place le socle du système pour éviter tout déplacement ultérieur (reportez-vous à Verrouillage du socle du système (▶ page 119)).

5.8.8 Diriger le câble de signal dans les serre-câbles

Il est possible d'utiliser les serre-câbles situés à l'arrière du système pour guider les câbles de signal et d'alimentation de manière organisée. La pompe et le compartiment à colonnes en option sont chacun munis d'un serre-câbles de chaque côté, tandis que le passeur d'échantillon possède deux serre-câbles de chaque côté. L'image indique l'emplacement des serre-câbles.



Illustration 67: Serre-câbles dans le système

5.9 Connexion d'un spectromètre de masse

Éléments nécessaires

Câble de signalement numérique d'entrée et de sortie



Illustration 68: Câble de signalement numérique d'entrée et de sortie

Broche	Description
1	Connecteur à 6 broches, pour la connexion au port Dig I/O du système
2	6 fils, pour la connexion à un appareil externe, par exemple, un spectromètre de masse

Procédez comme suit

- 1. Sur le passeur d'échantillon, branchez le connecteur à 6 broches du câble sur le port étiqueté **Dig I/O 2**.
- Branchez les fils pour les signaux de sortie (fil blanc et fil brun) sur les connecteurs appropriés du spectromètre de masse. Cette connexion permet d'obtenir automatiquement le signal de réponse d'injection.

5.10 Verrouillage du socle du système

Le socle du système a été mis en mode de déplacement après le déballage du système. Les outils de verrouillage se trouvent donc déjà en place dans les orifices du socle du système.

- 1. Placez le système dans la position d'utilisation.
- 2. Verrouillez le socle du système en faisant tourner les outils de verrouillage comme illustré ci-dessous.



Illustration 69: Verrouillage du socle du système (vue de haut)

3. Retirez les outils de verrouillage et rangez-les dans un endroit sûr, le tiroir du socle du système par exemple.

5.11 Mise en place des raccordements fluidiques

Le système est expédié avec la plupart des raccordements fluidiques installés pour un flux de travail d'injection directe. Suivez les instructions de la présente section pour compléter la configuration de tous les raccordements fluidiques.

NOTE Pour passer à un autre flux de travail, complétez l'installation tel que décrit dans le présent manuel, et exécutez le script : **A06 – Modifier le circuit fluidique / flux de tâches**. Pour en savoir plus sur d'autres flux de travail, consultez le *Guide de l'utilisateur du système Vanquish Neo*.

5.11.1 Passage des capillaires et des tubes dans le système

Les raccordements fluidiques entre les modules du système Vanquish transitent par la gaine de tubulure située à l'intérieur du système ou par les trous de guidage ou les attaches pour capillaires.

Gaine munie de guides tubulure

L'intérieur des modules empilables est doté d'une gaine de tubulure, sur le côté droit, qui permet de passer certains tubes et conduites depuis le module supérieur jusqu'au module inférieur, dans le système empilé Vanquish Neo. Cette gaine de tubulure est munie de quatre guides de tubulure.

Enfoncez les tubes ou les conduites dans les guides de tubulure.



Illustration 70: Gaine de tubulure munie de guides de tubulure (à gauche : vue depuis l'intérieur ; à droite : vue depuis le dessus)

Supports de tubulure

Des supports permettent de maintenir la tubulure en place. Glissez le côté du support sur la conduite d'évacuation.



Illustration 71: Support de tubulure (à gauche), support de tubulure installé (à droite)

Trous de guidage et attaches pour capillaire

Les trous de guidage et attaches pour capillaire sont fournis à des positions spécifiques sur les modules du système. Dans le système Vanquish Neo, dirigez les raccordements fluidiques d'un module au module suivant, en utilisant le trou de guidage ou l'attache pour capillaire correspondant, conformément aux instructions du manuel.

5.11.2 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.

5.11.2.1 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.

5.11.2.2 Connexion de capillaires nanoViper

La présente section décrit comment connecter les capillaires nanoViper[™] (ci-après dénommés capillaires Viper[™]).



Illustration 72: Connecteur Viper

Les connecteurs Viper sont installés sans molette. Si nécessaire, vous pouvez utiliser les molettes du kit d'expédition.

AVIS

- Serrez ou desserrez les capillaires Viper *uniquement* avec vos doigts et avec l'outil dynamométrique spécifique. N'utilisez pas d'outils autres que ceux fournis par Thermo Fisher Scientific à cette fin. Utilisez cet outil uniquement pour les raccordements pour lesquels il est prévu.
- 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.

Outils nécessaires

Tournevis dynamométrique comprenant :

- Base
- Adaptateur d'extension
- Mèche Viper

Procédez comme suit

AVIS—Endommagement du raccordement nanoViper

Attrapez le tournevis dynamométrique par sa base et maintenez-le droit. Si vous penchez le tournevis, une tension trop forte sera appliquée sur le raccordement nanoViper, ce qui pourrait endommager le raccordement ou le tournevis.

- 1. Insérez le capillaire Viper dans l'orifice de raccordement.
- 2. Serrez le raccord avec vos doigts.
- 3. Utilisez le tournevis pour serrer davantage le raccord. Pour accéder aux capillaires difficiles d'accès, utilisez l'adaptateur d'extension.



Illustration 73: Tournevis dynamométrique

N°	Description
1	Mèche Viper
2	Adaptateur d'extension
3	Base

- 4. Serrez les raccordements jusqu'à la limite de couple, signalée par un clic.
- 5. Vérifiez que le raccord ne fuit pas. S'il fuit, réparez la fuite.

Résolution de fuites

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

5.11.3 Raccordement des conduites d'évacuation

5.11.3.1 Vue d'ensemble des raccordements d'évacuation



AVERTISSEMENT—Fuite de substances dangereuses

Si les raccords des conduites d'évacuation et les bacs d'évacuation ne sont pas correctement installés, des substances dangereuses peuvent s'en échapper. L'exposition à ces substances peut présenter des risques pour la santé et la sécurité.

Vérifiez la bonne installation des raccordements et des bacs en testant le dispositif d'évacuation (reportez-vous à Vérification du dispositif d'évacuation () page 128)).

En cas de fuite dans le système, le liquide est évacué par l'orifice du dispositif d'évacuation du système situé sur le socle du système Vanquish.

Évacuation des liquides

Les liquides suivants s'écoulent par l'orifice d'évacuation situé en bas à droite sur le bac de fuite pour ensuite être récupérés par le dispositif d'évacuation :

- les liquides de fuite qui s'accumulent dans les bacs de fuite ;
- les déchets liquides provenant des dispositifs de rinçage de l'aiguille et la pompe d'évacuation dans le passeur d'échantillon;
- le déchet liquide provenant du débitmètre.



Illustration 74: Vue d'ensemble de la conduite d'évacuation

N°	Description
1	Conduite d'évacuation de rinçage de l'aiguille
2	Tubes d'évacuation de la pompe d'évacuation
3	Conduite d'évacuation du débitmètre (Purge Out)

Pour acheminer les liquides vers l'évacuation, reportez-vous à la section Raccordement du tube d'évacuation à l'orifice d'évacuation du système (▶ page 126).

Pour tester le dispositif d'évacuation, quand la conduite d'évacuation est installée, reportez-vous à la section Vérification du dispositif d'évacuation (> page 128).

NOTE Si vous devez raccourcir la longueur du tube, utilisez un coupetube. 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.

5.11.3.2 Raccordement du tube d'évacuation à l'orifice d'évacuation du système

Éléments nécessaires

- Tube d'évacuation du kit d'expédition
- Raccord coudé du kit d'expédition
- Bac d'évacuation approprié

Procédez comme suit

1. Insérez aussi loin que possible le raccord coudé dans l'orifice d'évacuation du système pour obtenir un raccord étanche.



Illustration 75: Tube d'évacuation du système et raccord coudé connectés

N°	Description
1	Pied d'évacuation du système
2	Sortie d'évacuation du détecteur
2a	Conduite d'évacuation du détecteur
3	Raccord coudé
3a	Conduite d'évacuation du système

- 2. Insérez manuellement aussi loin que possible le tube d'évacuation dans le raccord coudé pour obtenir un raccord étanche.
- 3. Dirigez l'extrémité libre du tube d'évacuation vers un bac d'évacuation.

4. Uniquement si le tube d'évacuation est dirigé vers l'arrière du système : introduisez le tube d'évacuation dans les clips et assurezvous qu'aucun siphon n'a été formé.



Illustration 76: Insertion du tube d'évacuation dans les clips

5. Assurez-vous que les déchets s'évacuent correctement (reportezvous à la section suivante).

Vérification de l'évacuation des déchets

- 1. Placez le bac d'évacuation plus bas que le système empilé. Assurezvous que la distance est suffisante pour empêcher les vapeurs de solvant de retourner dans le système.
- 2. Assurez-vous que le tube d'évacuation n'est ni plié ni pincé.
- Pour empêcher le liquide de refluer dans le système, l'ensemble du tube d'évacuation doit se trouver plus bas que l'orifice d'évacuation du système. Veillez à éviter la formation de siphons.
- Pour permettre au liquide évacué de s'écouler correctement et prévenir toute accumulation de liquide dans le socle du système, l'extrémité du tube d'évacuation ne doit pas baigner dans le liquide évacué (reportez-vous à la figure ci-dessous) :
 - a) Si nécessaire, raccourcissez le tube d'évacuation.
 - b) Surveillez le niveau de liquide dans le bac d'évacuation et videzle au besoin.



Illustration 77: Extrémité libre du tube d'évacuation dans le bac d'évacuation

5.11.3.3 Vérification du dispositif d'évacuation

Vérifiez que le dispositif d'évacuation fonctionne correctement :

- 1. Versez de l'eau ou de l'isopropanol dans l'évacuation du bac à solvants.
- 2. Vérifiez que le liquide s'évacue du système correctement par l'orifice d'évacuation et s'écoule dans le bac d'évacuation.
- 3. En cas de fuite, trouvez et corrigez le problème.

5.12 Remplissage des réservoirs de solvant

Avant d'utiliser le système, remplissez les réservoirs de solvant.

NOTE Pour éviter la formation de bulles d'air dans les conduites de liquide, placez toujours les réservoirs dans le bac à solvant sur le dessus du système empilé.

Préparatifs

- 1. Rincez soigneusement les réservoirs de solvant et toutes les pièces du filtre de conduite de solvant. Utilisez un solvant ultrapur pour le rinçage.
- 2. Montez le filtre de conduite de solvant en revêtant des gants propres et appropriés :
 - a) Insérez l'adaptateur de conduite de solvant dans le porte-filtre (partie supérieure).
 - b) Tournez le porte-filtre (partie supérieure) avec adaptateur tête en bas et pressez-le sur une surface appropriée pour que l'adaptateur soit fermement attaché sur le porte-filtre.
 - c) Placez le fritté dans une position de niveau dans le porte-filtre (partie inférieure).
 - d) Vissez le haut du filtre fermement sur le bas du filtre.



Illustration 78: Filtre de conduite de solvant

N°	Description
1	Adaptateur de conduite de solvant
2	Porte-filtre, partie supérieure
3	Fritté de filtre
4	Porte-filtre, partie inférieure

Procédez comme suit



Illustration 79: Préparation de la conduite de solvant et du bouchon de réservoir

N°	Description
1	Conduite de solvant
2	Guide de maintien
3	Bouchon de réservoir
4	Capuchons protecteurs
5	Filtre de conduite de solvant avec adaptateur de conduite de solvant

1. Faites passer chaque conduite de solvant par un guide de maintien.

- Faites passer la conduite de solvant par un trou dans un bouchon pour réservoir de solvant. Le guide de maintien permet d'immobiliser la conduite de solvant dans le réservoir.
- 3. Obturez les trous restants des bouchons de réservoir au moyen de capuchons protecteurs.
- 4. Glissez un filtre de conduite de solvant sur chaque conduite de solvant.
- 5. Remplissez les réservoirs de solvant avec du solvant.
- 6. Serrez le bouchon du réservoir manuellement. Enfoncez le guide de maintien dans le trou du bouchon du réservoir afin d'immobiliser la tubulure dans le bouchon.
- 7. Placez les réservoirs de solvant dans le bac à solvants.
- 8. Positionnez les conduites de solvant directement dans les guides de tubulure.
- Contrôlez les conduites de solvant sur l'ensemble du trajet d'écoulement. Assurez-vous que les conduites ne sont, à aucun endroit du trajet d'écoulement, pliées, pincées ou serrées.

Consultez également

Informations sur les solvants et les additifs (> page 144)

5.13 Remplissage des réservoirs de rinçage de l'aiguille

Avant d'utiliser le système, vous devez remplir les réservoirs de rinçage de l'aiguille et purger les dispositifs de rinçage de l'aiguille.

NOTE Pour éviter la formation de bulles d'air dans les conduites de liquide, placez toujours les réservoirs dans le bac à solvant sur le dessus du système empilé.

Outils nécessaires

Coupe-tube

Préparatifs

Rincez soigneusement les réservoirs de rinçage. Utilisez un solvant ultrapur pour le rinçage.

Liquides de rinçage de l'aiguille

Sélectionnez des liquides de rinçage de l'aiguille adaptés à votre application et qui éliminent suffisamment les échantillons résiduels de l'aiguille. Il y a un réservoir de liquide de rinçage fort et un autre de liquide de rinçage faible. Généralement, le liquide de rinçage de l'aiguille est le solvant B pour votre application ou un solvant plus fort, tandis que le liquide de rinçage faible est proche du solvant de début de votre gradient.

Type de liquide de rinçage de l'aiguille	Description (pour les applications en phases inversées)	Tubes
Fort	Liquide de rinçage riche en matière organique, par exemple, 0,1 % d'acide formique dans de l'acétonitrile	Conduite de rinçage pour purge et arrêt de courte durée, étiquetée S (rouge) : connectée à Strong l'orifice d'admission du dégazeur situé sur la pompe.
		Conduite de rinçage pour le rinçage de l'intérieur et de l'extérieur de l'aiguille, étiquetée SWP (rouge) : connectée au port d'entrée supérieur de dégazeur sur la pompe de rinçage de l'aiguille.
Faible	Liquide de rinçage de l'aiguille faible en matière organique, par exemple, 0,1 % d'acide formique dans de l'eau	Conduite de rinçage pour le rinçage de l'intérieur de l'aiguille et pour purge, étiquetée W (verte) : connectée à Weak l'orifice d'admission du dégazeur situé sur la pompe.
		Conduite de rinçage pour le rinçage de l'extérieur de l'aiguille, étiquetée WWP (verte) : connectée au port d'entrée inférieur de dégazeur sur la pompe de rinçage de l'aiguille.

Procédez comme suit pour chaque réservoir



Illustration 80: Préparation de la conduite de rinçage et du bouchon de réservoir

N°	Description
1	Conduite de rinçage
2	Guide de maintien
3	Bouchon de réservoir
4	Capuchons protecteurs
5	Filtre de conduite de solvant avec adaptateur de conduite de solvant

- 1. Faites passer chaque conduite de rinçage de l'aiguille par un guide de maintien.
- 2. Faites passer chaque conduite de rinçage de l'aiguille par un trou du bouchon du réservoir de rinçage. Le guide de maintien permet d'immobiliser la conduite de solvant dans le réservoir.
- 3. Obturez les trous restants des bouchons de réservoir au moyen de capuchons protecteurs.
- 4. Glissez un porte-filtre avec adaptateur sur chaque conduite de rinçage.
- 5. Remplissez le réservoir de rinçage avec du liquide de rinçage.
- 6. Serrez le bouchon du réservoir manuellement. Enfoncez le guide de maintien dans le trou du bouchon du réservoir afin d'immobiliser la tubulure dans le bouchon.
- 7. Placez le réservoir de rinçage dans le bac à solvants.

- 8. Inspectez les conduites de rinçage sur l'ensemble du trajet d'écoulement :
 - Assurez-vous que les conduites de rinçage ne sont, à aucun endroit du trajet d'écoulement, pliées (plicaturées), pincées ou serrées.
 - b) Si vous devez raccourcir la longueur du tube, utilisez un coupetube. Veillez à ce que la coupe soit à angle droit par rapport à l'axe de la conduite.

Purge des dispositifs de rinçage de l'aiguille

Les dispositifs de rinçage de l'aiguille sont purgés quand vous exécutez le script **A02 – Démarrer automatiquement** après avoir mis le système en marche. Reportez-vous à la section Préparation du système pour l'utilisation initiale () page 139).

5.14 Remplissage du réservoir de liquide de rinçage des joints

Avant d'utiliser le système, remplissez le réservoir de liquide de rinçage des joints.

NOTE Pour éviter la formation de bulles d'air dans les conduites de liquide, placez toujours les réservoirs dans le bac à solvant sur le dessus du système empilé.

Outils nécessaires

Coupe-tube

Préparatifs

Rincez soigneusement le réservoir de liquide de rinçage des joints. Utilisez un solvant ultrapur pour le rinçage.

Liquide de rinçage des joints

Utilisez de l'isopropanol à 75 % dans de l'eau avec de l'acide formique à 0,1 % (qualité HPLC ; préparation en volume, par exemple, 75 mL d'isopropanol + 25 mL d'eau + 0,1 mL d'acide formique).

Procédez comme suit





N°	Description
1	Conduite de rinçage
2	Guide de maintien
3	Bouchon de réservoir
4	Capuchons protecteurs

- 1. Faites passer la conduite de rinçage par le guide de maintien.
- Faites passer la conduite de rinçage par un trou du bouchon du réservoir de rinçage. Le guide de maintien permet d'immobiliser la tubulure dans le réservoir.
- 3. Obturez les trous restants du bouchon de réservoir au moyen de capuchons protecteurs.
- 4. Remplissez le réservoir de rinçage avec du liquide de rinçage.
- 5. Serrez le bouchon du réservoir manuellement. Enfoncez le guide de maintien dans le trou du bouchon du réservoir afin d'immobiliser la tubulure dans le bouchon.
- 6. Placez le réservoir de rinçage dans le bac à solvants.
- 7. Inspectez les conduites de rinçage sur l'ensemble du trajet d'écoulement :
 - Assurez-vous que les conduites de rinçage ne sont, à aucun endroit du trajet d'écoulement, pliées (plicaturées), pincées ou serrées.
 - b) Si vous devez raccourcir la longueur du tube, utilisez un coupetube. Veillez à ce que la coupe soit à angle droit par rapport à l'axe de la conduite.

5.15 Première mise en marche du système

Si la configuration de votre système comprend un compartiment à colonnes, installez-le avant de mettre le système en marche.

Boutons d'alimentation et interrupteurs d'alimentation principaux

L'image illustre les dispositifs d'alimentation du système.



Illustration 82: Boutons d'alimentation et interrupteurs d'alimentation principaux

N°	Description
1	Bouton d'alimentation du système
2	Bouton d'alimentation du contrôleur du système
3	Interrupteurs d'alimentation principaux

Préparatifs

1. Vérifiez que le verrou d'expédition du passeur d'échantillon est correctement desserré.

AVIS—Dommages au système

Avant de mettre le système Vanquish sous tension, vérifiez que le verrou d'expédition a été correctement desserré afin d'éviter d'endommager le système.

 Vérifiez que le logiciel de chromatographie est installé sur l'ordinateur du système de données. Quand l'appareil est mis sous tension, les pilotes USB nécessaires sont recherchés automatiquement et le système d'exploitation Windows™ peut alors détecter les modules du système.

Procédez comme suit

- Mettez chaque module du système sous tension (pompe, passeur d'échantillon, compartiment à colonnes en option) au moyen de son interrupteur d'alimentation principal. Mettez le passeur d'échantillon sous tension avant de démarrer le débit de la pompe et avant que la pression ne monte. La pompe démarre et contrôle le système de rinçage des joints. Chaque fois que vous mettez la pompe sous tension, elle commence son fonctionnement avec un cycle de rinçage des joints.
- Mettez le contrôleur du système sous tension au moyen de son interrupteur d'alimentation principal.
 Il suffit d'appuyer sur l'interrupteur d'alimentation principal pour mettre le contrôleur du système sous tension et il n'est pas nécessaire d'utiliser le bouton-poussoir d'alimentation. Après le démarrage du contrôleur du système, la lumière LED d'état s'allume.
- 3. Mettez le système sous tension au moyen du bouton d'alimentation électrique du système.

5.16 Intégration dans un appareil LAN

Pour activer la communication entre le système et un appareil LAN, y compris l'ordinateur des données chromatographiques du système et un autre dispositif, comme un spectromètre de masse, vous devez l'intégrer dans un appareil LAN. La Vanquish User Interface propose deux différentes possibilités de connectivité :

- au moyen d'une adresse IP personnalisée Requis, par exemple, si un dispositif dans l'appareil LAN possède une adresse IP personnalisée.
- au moyen d'un protocole de configuration dynamique des hôtes (DHCP)

Prérequis

- Vous avez branché le contrôleur du système sur l'ordinateur.
- Vous avez configuré un appareil LAN sur un ordinateur sur lequel est installé le système de gestion de données chromatographiques.
- Pour la possibilité d'adresse IP personnalisée seulement : Vous avez obtenu une plage d'adresses IP et le masque de réseau de l'appareil LAN.

Procédez comme suit

- 1. Sur le dispositif d'affichage, ouvrez le volet paramètres.
- 2. Sélectionnez la boîte **Connectivité**.
- 3. Sélectionnez l'une des possibilités de connectivité.
- 4. Si vous avez sélectionné Personnalisé :
 - a) Saisissez le masque de réseau de l'appareil LAN.
 - b) Saisissez une adresse IP unique du contrôleur du système dans votre appareil LAN.
- 5. Sélectionnez Appliquer.

Consultez également

Connexion d'un câble LAN (> page 115)

5.17 Préparation du système pour l'utilisation initiale

AVIS—Les particules et les bulles d'air peuvent nuire au fonctionnement du système

Certains composants du système sont remplis d'isopropanol quand ils sont entreposés ou quand ils sont expédiés du site de fabrication.

- Pour éliminer les bulles d'air (et l'isopropanol) avant d'utiliser le système pour la première fois, exécutez le script : A02 – Démarrer automatiquement.
- Pour éliminer l'isopropanol, utilisez des solvants miscibles à l'isopropanol. S'ils ne le sont pas, utilisez un solvant intermédiaire approprié.

À quel moment exécuter cette opération ?

Avant la toute première utilisation du système

Liquide requis	Consommation pendant l'exécution du script : A02 – Démarrer automatiquement	Consommation pendant l'analyse	Exigences de liquide
Solvant	Max. 10 mL par réservoir de solvant	En fonction du débit	Lorsque vous utilisez le système pour la première fois, employez des solvants miscibles à l'isopropanol. S'ils ne le sont pas, utilisez un solvant intermédiaire approprié. Pour tous les autres cas, reportez-vous à la section Informations sur les solvants et les additifs () page 144).
Liquide de rinçage des joints	Max. 10 mL	Environ 250 mL tous les 14 jours	Reportez-vous à la section Remplissage du réservoir de liquide de rinçage des joints () page 134).

Quantités requises de solvant et de liquide de rinçage des joints

Consommation pendant le script : A02 – Démarrer automatiquement	Consommatic l'analyse	on pendant	Exigences de liquide	
Max. 10 mL	Consommation par réservoir de rinçage de l'aiguille et par injection ¹ :		Reportez-vous à la section Remplissage des réservoirs de rinçage de l'aiguille	
	SWP	500 μL	(▶ page 131).	
	WWP	400 μL	-	
	W	250 μL		
¹ : Les valeurs s'appliquent à : boucle d'échantillonnage de 25 µL, injection directe				

Quantités requises de liquide de rinçage de l'aiguille

¹ : Les valeurs s'appliquent à : boucle d'échantillonnage de 25 μL, injection directe (boucle en bypass) configurée dans le système de gestion de données chromatographiques, réglages de rinçage par défaut du système de gestion de données chromatographiques.

Remarque : la consommation dépend de la taille de la boucle d'échantillonnage, la taille de la colonne, la durée du rinçage, le flux du travail et la vitesse de rinçage. Si vous utilisez d'autres réglages, les valeurs pourraient être significativement plus élevées.

Procédez comme suit

- 1. Vérifiez le niveau de liquide de tous les réservoirs, surtout celui des réservoirs de rinçage de l'aiguille. Assurez-vous que les quantités de liquide sont suffisantes pour les points suivants :
 - le script A02 Démarrer automatiquement (reportez-vous au tableau ci-dessus);
 - la première analyse (reportez-vous au tableau ci-dessus et ajoutez les quantités selon le débit de l'application).
- 2. Fermez les portes de tous les modules du système Vanquish Neo, si ce n'est pas déjà fait.
- Pour configurer le solvant A/B, exécutez le script : A01 Définir le type de solvant de la pompe.
- Exécutez le script avec ou sans diagnostic (test de fuite) : A02 Démarrer automatiquement.
 Le script rince le trajet d'écoulement du système jusqu'au capillaire d'entrée de la colonne, teste la contre-pression du système et effectue un test de fuite, le cas échéant.
- Pour configurer un type de colonne, exécutez le script : A03 Définir le type de colonne de séparation.
- 6. Pour configurer les propriétés de la colonne, exécutez le script : A04
 Définir les caractéristiques de la colonne de séparation.
- Pour préparer la colonne, exécutez le script : B06 Conditionnement de la / des colonne(s).

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 Consignes de sécurité pendant l'utilisation

Lorsque vous utilisez le système, 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 26).



AVERTISSEMENT—Pièces mobiles

Les pièces internes du passeur d'échantillon se déplacent lorsque le passeur d'échantillon se prépare et procède aux analyses d'échantillon. Ces pièces mobiles peuvent présenter un risque de pincement et entraîner des blessures corporelles.

- Maintenez les portes avant du passeur d'échantillon fermées lorsque le passeur d'échantillon se prépare et procède aux analyses d'échantillon.
- Durant ces phases, la barre à LED située sur le passeur d'échantillon s'illumine en bleu.



AVERTISSEMENT—Vapeurs inflammables et dangereuses provenant de déversements

Des vapeurs inflammables ou dangereuses provenant de déversements d'échantillons peuvent s'accumuler à l'intérieur du passeur d'échantillon. Cette accumulation peut présenter des risques pour la santé et la sécurité.

- Assurez-vous que les portoirs d'échantillons et plaques à puits sont correctement placés dans les segments.
- Si vous utilisez un portoir d'échantillons, remplissez-le avec les flacons avant de le placer dans le carrousel.
- En cas de déversement à l'intérieur du passeur d'échantillon, mettez-le hors tension. Nettoyez le déversement et laissez la porte du passeur d'échantillon ouverte. Patientez le temps nécessaire au séchage du déversement et à la dispersion des vapeurs avant de reprendre l'utilisation du passeur d'échantillon.



ATTENTION—Surfaces chaudes

Les surfaces internes du système 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.

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.
- Vérifiez toujours que le passeur d'échantillon est allumé avant de mettre la pompe sous pression et d'augmenter la pression du système. Si le passeur d'échantillon est hors tension (par exemple, à la suite d'une coupure de courant), arrêtez la pompe et patientez jusqu'à ce que la pression redescende à zéro avant de rallumer le passeur d'échantillon ou d'autres modules.

6.2 Informations sur les solvants et les additifs

6.2.1 Compatibilité générale

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

- Utilisez uniquement des solvants et additifs compatibles avec tous les composants du trajet d'écoulement. Pour obtenir des informations sur les matériaux utilisés dans le trajet d'écoulement analytique, reportez-vous à la section Caractéristiques de performance (> page 401) de chaque module du système.
- L'insertion de bouteille est composée de polypropylène expansé à cellules fermées, qui est compatible avec la plupart des solvants HPLC communs.
- Le méthanol anhydre peut mener à des défaillances de stress dans les surfaces en titane, surtout quand de l'acide formique est ajouté au mélange. Thermo Fisher Scientific recommande d'ajouter 3 % d'eau pour éviter que cela ne survienne.

Compatibilité avec le joint de piston

- La pompe est livrée avec des joints de piston en PE-UHMW.
 L'utilisation de tétrahydrofurane, de cétones ou d'hydroxyde d'ammonium comme solvants peut provoquer une détérioration des joints.
- 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.

6.2.2 Plages pH autorisées

Plage pH autorisée pour le système :

Plage pH autorisée	Remarque
2-10	Valeurs de pH de 2 : La durée d'application doit être aussi brève que possible. Rincez abondamment le système après ces applications.
6.2.3 Concentrations autorisées

Concentration	Valeur
Concentration tampon	s.o.
Concentration en chlorure	Inférieure ou égale à 1 mol/l Avec une concentration élevée en chlorure, la durée d'application doit être aussi brève que possible. Rincez abondamment le système après ces applications.

Concentrations autorisées pour le système :

6.2.4 Utilisation de solvants et d'additifs

Les particules qui pénètrent dans le système chromatographique peuvent obstruer les capillaires et les vannes, augmenter l'usure et endommager la colonne ou le système. Des algues et autres microorganismes peuvent se développer et se déposer dans le système chromatographique, colmatant ainsi les filtres des conduites de solvant, en particulier lors de l'utilisation de solvants aqueux. L'obstruction de capillaires ou de filtres peut provoquer une augmentation ou des fluctuations de la pression du système.

Pour un fonctionnement optimal du système, respectez les consignes suivantes.

Qualité de solvant et filtration

- Utilisez des solvants et additifs ultrapurs de qualité élevée (filtrés) conformément aux exigences de l'application, par exemple de qualité UHPLC ou LC/MS. Les solvants ultrapurs filtrés sont généralement étiquetés de cette manière par le fournisseur.
- Utilisez de l'eau de qualité élevée, par exemple, de niveau UHPLC ou LC/MS (filtration à 0,2 μm).
- Lorsque vous utilisez de l'eau provenant de systèmes de purification, elle peut être contaminée par des polymères si le système n'est pas correctement entretenu.
- Lorsque la préparation de solutions salines ou de tampons est terminée, filtrez-les en utilisant une membrane (0,2 μm) afin de retirer toute particule et de diminuer la croissance microbienne.
- Utilisez les filtres de conduite de solvant appropriés recommandés par Thermo Fisher Scientific. Vérifiez régulièrement la perméabilité des frittés de filtres et remplacez-les le cas échéant.

- Utilisez des solvants frais correctement dégazés. Remplacez les solvants au moins une fois par semaine. Remplacez toujours tout le solvant dans le réservoir. Ne rajoutez pas simplement du solvant, car cela pourrait causer des changements de composition du solvant ou la formation de composants non souhaités dans les réservoirs.
- Pour prolonger la durée de vie des joints, ne recyclez pas le solvant et ne le faites pas circuler dans le système de nouveau.
- Notez bien les propriétés propres à chaque solvant, comme la viscosité, la miscibilité, le point d'ébullition ou l'absorption des UV, ainsi que la solubilité des sels de toutes les proportions pendant le gradient.

Réservoirs de solvant

- Avant de remplir un réservoir de solvant, rincez-le soigneusement à l'aide d'un solvant ultrapur.
- Pour réduire le développement des algues, utilisez des flacons en verre ambré. Utilisez des additifs appropriés, par exemple l'acide formique, pour les solvants aqueux.
- Évitez d'utiliser des détergents quand vous nettoyez les flacons.
 Rincez les flacons utilisés pour les applications LC/MS (y compris les cylindres gradués) avec des solvants de qualité LC/MS avant de les utiliser, et étiquetez-les et entreposez-les séparément.
- N'utilisez pas du méthanol provenant de réservoirs en aluminium.

Rinçage après utilisation

- Rincez le système pour éliminer les tampons et solutions susceptibles de former des peroxydes.
- Ne laissez jamais de tampons, de solutions salines ou de solvants agressifs dans le système pendant une période prolongée sans les faire circuler.
- Avant de passer d'une solution saline ou tampon à un solvant organique, rincez soigneusement le système avec de l'eau de qualité UHPLC ou LC/MS.
- Avant de passer à un autre type de solvant, assurez-vous que le nouveau solvant est miscible avec le solvant précédent. S'il ne l'est pas, une floculation peut se produire.
- Basculez entre deux solvants immiscibles en passant par un solvant intermédiaire, par exemple, de l'isopropanol. Étape par étape, remplacez le premier solvant par le solvant intermédiaire, puis remplacez, étape par étape, le solvant intermédiaire par le nouveau solvant.

6.3 Consignes générales pour l'utilisation du système

Respectez l'information et les consignes suivantes lors de l'utilisation du système.

Pour une performance optimale

- Rincez l'intérieur et l'extérieur de l'aiguille pour réduire le transfert entre échantillons. Le rinçage de l'aiguille permet de s'assurer que tout échantillon résiduel est correctement éliminé de l'aiguille.
- Gardez la porte du compartiment à colonnes fermées pendant cette opération.
- Évitez d'ouvrir les vannes à sec, c'est-à-dire sans débit de pompe. Cela peut détériorer la vanne.
- N'allumez pas la pompe lorsque le système de rinçage des joints est sec.
- Respectez les consignes générales et les recommandations sur l'utilisation de solvants et d'additifs dans le système. Reportez-vous à la section Informations sur les solvants et les additifs (> page 144).

Notes concernant le traitement des échantillons

Notez bien les consignes suivantes quand vous utilisez le passeur d'échantillon :

NOTE—Garder les portes du passeur d'échantillon fermées pendant le traitement des échantillons

- Si vous ouvrez les portes pendant le déplacement de l'aiguille et du carrousel, le passeur d'échantillon interrompt le déplacement en cours du bras de l'aiguille. L'aiguille ne perce pas le récipient d'échantillon et ne prélève pas l'échantillon. Un signal sonore est émis pour vous avertir de l'ouverture des portes.
- Chaque fois que les portes sont fermées, le passeur d'échantillon inventorie les portoirs d'échantillons et les récipients d'échantillonnage présents dans le compartiment à échantillons.

Consultez également

Optimisation des performances (> page 169)

6.4 Mise en marche du système

- Mettez le contrôleur du système en marche en appuyant sur le bouton pressoir supérieur sur le socle du système. Après le démarrage du contrôleur du système, la lumière LED d'état s'allume.
- 2. Mettez le système sous tension au moyen du bouton d'alimentation électrique inférieur sur le socle du système.

6.5 Éléments de commande

Le système est conçu pour être piloté principalement depuis un ordinateur équipé du logiciel de chromatographie et de Vanquish User Interface.

De plus, les éléments suivants sont également disponibles dans le système :

- Panneaux de commande :
 Les boutons du panneau de commande vous permettent d'exécuter certaines fonctions directement depuis l'appareil.
- Voyants d'état : Permettent un contrôle visuel rapide de l'état de fonctionnement de l'appareil. Le système est muni des voyants d'état suivants :
 - Barre à LED de voyants d'état située sur le devant de chaque appareil.
 - **STATUS** LED sur les panneaux de commande.
 - LED d'état du contrôleur du système

6.5.1 Panneaux de commande

Les éléments suivants sont disponibles :

STATUS LED

La LED **STATUS** permet un contrôle visuel rapide de l'état opérationnel de l'appareil. Lorsque les portes sont fermées, la barre à LED située sur la façade avant indique l'état opérationnel.

• Boutons

Les boutons vous permettent d'exécuter certaines fonctions directement depuis l'appareil. Lorsque vous appuyez sur un bouton, un bref signal sonore confirme que la fonction s'exécute.

LED D'ÉTAT

La LED **STATUS** du panneau de commande 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. Vérifiez le message sur le volet notification. Pour connaître les mesures appropriées à prendre, reportez-vous à la section Dépannage (> page 353).

Pour obtenir des informations concernant les LED situées à côté d'un bouton sur le panneau de commande, reportez-vous au tableau Boutons du panneau de commande.

Boutons du panneau de commande

Les boutons ont les fonctions suivantes :

Passeur d'échantillon et pompe

	Bouton	Fonction
	DOCK	Appuyer sur ce bouton permet de déplacer les pistons de la tête de pompe sélectionnée vers une position de retrait appropriée pour les procédures de maintenance, comme le remplacement de tête de pompe ou de piston.
		Quand la tête de pompe est dans la position de maintenance, la LED située à côté du bouton FLOW commence à clignoter en vert. En appuyant de nouveau sur le bouton, la tête de pompe retourne à la position d'utilisation. La LED arrête de clignoter en vert.
	FLOW	 Appuyer sur ce bouton permet de démarrer ou d'arrêter le débit, en tenant compte du débit sélectionné et de la composition de solvant. Les LED L et R à côté du bouton réfèrent aux têtes de pompe gauche ou droite : La LED est éteinte = le débit de la pompe est désactivé ou à zéro La LED est verte = la tête de pompe est en fonctionnement La LED clignote en vert = la tête de pompe est dans la position des procédures de maintenance
	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 dans les 10 minutes. Sinon, le signal sonore reprend. Si l'appareil détecte un problème différent, le signal sonore vous prévient de nouveau immédiatement.
	PURGE	Réservé pour utilisation ultérieure. Pour de l'information sur la purge, reportez-vous à la section Purge (vue d'ensemble) (> page 65).
	ROTATE	Appuyer sur ce bouton permet de faire tourner le carrousel dans le sens antihoraire jusqu'à la position de chargement suivante.
	SELECT	 Passeur d'échantillon Réservé pour utilisation ultérieure. Pompe Appuyer sur ce bouton permet de sélectionner la tête de pompe pour laquelle vous pouvez lancer la fonction DOCK en utilisant le panneau de commande : La LED L est verte = la tête de pompe gauche est sélectionnée La LED R est verte = la tête de pompe dreite est sélectionnée
		• La LED K est verte = la tete de pompe droite est selectionnée

Bouton	Fonction
SERVICE	Le bouton a les fonctions suivantes :
	 Appuyer une fois sur ce bouton lève l'aiguille du siège de l'aiguille et règle la vanne d'injection en position Bypass. Cette opération est nécessaire par exemple pour serrer la vis de verrouillage d'expédition et fixer l'unité d'aiguille en place.
	 Appuyer sur ce bouton une deuxième fois initialise l'aiguille et elle se déplace au-dessus de l'orifice de rinçage.
	 Appuyer une troisième fois sur ce bouton fait redescendre l'aiguille dans son siège et initialise l'aiguille. Remarque : La vis de verrouillage d'expédition doit être desserrée.
VALVE	Appuyer sur ce bouton permet de basculer la vanne d'injection en position Bypass . La LED B est verte lorsque la vanne d'injection est dans la position Bypass .
WASH	Appuyer sur ce bouton lance un cycle de rinçage extérieur de l'aiguille au cours duquel la surface externe de l'aiguille est rincée dans l'orifice de rinçage. Après le cycle de rinçage, l'aiguille retourne dans son siège.

Compartiment à colonnes

Bouton	Fonction
MUTE ALARM	Un signal sonore vous prévient lorsque l'appareil détecte un problème. Pour désactiver le signal sonore de l'alarme en cours, appuyez sur ce bouton. Éliminez la source de l'alarme dans les 10 minutes. Sinon, le signal sonore reprend. Si l'appareil détecte un problème différent, le signal sonore vous prévient de nouveau immédiatement.
TEMP CTRL	 Appuyer sur ce bouton allume ou arrête la régulation de la température : La LED à côté du bouton est verte = la régulation de la température est activée
	 La LED a cote du bouton est éteinte = la regulation de la température est désactivée
VALVE LEFT	Appuyer sur ce bouton permet d'activer la vanne de commutation pour colonne gauche (si installée).
	Garder ce bouton enfoncé active la fonction de vannes de capture si une vanne VICI est installée dans la position de vanne gauche.
	Pour les vannes à 2 positions et 6 ports, les LED situées à côté des boutons indiquent la position de vanne actuelle :
	• La LED du haut est verte = la vanne est dans la position 1
	• La LED du bas est verte = la vanne est dans la position 2
	 Les LED sont éteintes = la vanne bascule

Bouton	Fonction
VALVE RIGHT	Appuyer sur ce bouton permet d'activer la vanne de commutation pour colonne droite (si installée).
	Garder ce bouton enfoncé active la fonction de vannes de capture si une vanne VICI est installée dans la position de vanne droite.
	Pour les vannes à 2 positions et 6 ports, les LED situées à côté des boutons indiquent la position de vanne actuelle :
	 La LED du haut est verte = la vanne est dans la position 1
	• La LED du bas est verte = la vanne est dans la position 2
	 Les LED sont éteintes = la vanne bascule

6.5.2 Barres à LED

Les barres à LED sur le devant des modules offrent de l'information quand les modules sont fermées. Quand le système est connecté dans le système de gestion de données chromatographiques, il est possible que les barres à LED donnent moins d'information.

Passeur d'échantillon

Barre à LED	Description
Éteinte (sombre)	Le module est hors tension.
Lumière estompée	Les portes du module sont ouvertes.
Clignotement lent jaune	L'alimentation du module est activée ou le module est équilibré, mais le système n'est pas connecté au système de gestion de données chromatographiques.
Jaune	Le système est connecté au système de gestion de données chromatographiques, mais le module n'est pas équilibré.
Verte, clignotante	Le module s'équilibre. Si vous utilisez la thermostatisation du compartiment à échantillons, la température de thermostatisation n'est pas encore atteinte.
Verte	Le système est connecté au système de gestion de données chromatographiques et le module est équilibré, mais aucune donnée n'est acquise. Si vous utilisez la thermostatisation du compartiment à échantillons, la température de thermostatisation est atteinte.
Bleue, en cours	Le passeur d'échantillon prépare l'injection. Certaines pièces du passeur d'échantillon sont en mouvement.
Bleue	Le système est connecté au système de gestion de données chromatographiques et l'injection ou la séquence est en cours, y compris l'acquisition de données. Ou : Un test de diagnostic est en cours.
Rouge	Un problème ou une erreur est survenu. Vérifiez le message sur le volet notification. Pour connaître les mesures appropriées à prendre, reportez-vous à la section Dépannage () page 353).

Pompe

Barre à LED	Description
Éteinte (sombre)	Le module est hors tension.
Lumière estompée	Les portes du module sont ouvertes.
Clignotement lent jaune	L'alimentation du module est activée ou le module est équilibré, mais le système n'est pas connecté au système de gestion de données chromatographiques.

Barre à LED	Description
Jaune	Le système est connecté au système de gestion de données chromatographiques, mais le module n'est pas équilibré. Le débit de la pompe est arrêté.
Verte, clignotante	Un cycle de purge est en cours.
Verte	Le système est connecté au système de gestion de données chromatographiques et le module est équilibré, mais aucune donnée n'est acquise. Le débit de la pompe est allumé.
Bleue	Le système est connecté au système de gestion de données chromatographiques et l'injection ou la séquence est en cours, y compris l'acquisition de données. Ou : Un test de diagnostic est en cours.
Rouge	Un problème ou une erreur est survenu. Vérifiez le message sur le volet notification. Pour connaître les mesures appropriées à prendre, reportez-vous à la section Dépannage () page 353).

Compartiment à colonnes

Barre à LED	Description
Éteinte (sombre)	Le module est hors tension.
Lumière estompée	Les portes du module sont ouvertes.
Clignotement lent jaune	L'alimentation du module est activée ou le module est équilibré, mais le système n'est pas connecté au système de gestion de données chromatographiques.
Jaune	Le système est connecté au système de gestion de données chromatographiques, mais le module n'est pas équilibré.
Verte, clignotante	Le module s'équilibre.
Verte	Le système est connecté au système de gestion de données chromatographiques et le module est équilibré, mais aucune donnée n'est acquise.
Bleue	Le système est connecté au système de gestion de données chromatographiques et l'injection ou la séquence est en cours, y compris l'acquisition de données. Ou : Un test de diagnostic est en cours.
Rouge	Un problème ou une erreur est survenu. Vérifiez le message sur le volet notification. Pour connaître les mesures appropriées à prendre, reportez-vous à la section Dépannage (> page 353).

6.6 Vue d'ensemble de l'interface utilisateur Vanquish

6.6.1 Informations sur le contrôleur du système

Les informations sur le contrôleur du système s'affichent en haut à gauche.



Illustration 83: Informations du contrôleur du système

N°	Description
1	Heure ¹
2	Date ¹
3	Version du microprogramme du contrôleur du système
4	Numéro de série du contrôleur du système
¹ : Le format dépend de vos réglages de langue.	

6.6.2 Icônes d'état de l'état général du système

Les icônes suivantes représentent l'état général du système.

Icône	État général du système
	Ok
	Erreur Un ou plusieurs modules du système ont signalé une erreur et la barre à LED d'un ou de plusieurs modules est rouge.
	Pour obtenir des informations sur la résolution de l'état Erreur, reportez-vous à la section Dépannage du système (> page 379).

6.6.3 Icônes de la barre de navigation et autres icônes

La barre de navigation comprend les icônes suivantes.

lcône	Description
	Volet Accueil Donne accès aux paramètres fonctionnels de tous les modules du système connectés au contrôleur du système à l'aide de l'interface System Interlink. Pour obtenir des informations complémentaires, reportez-vous à la section Volet Accueil () page 157).
*	Volet script Offre des scripts automatisés. Pour obtenir des informations complémentaires, reportez-vous à la section Volet Script (▶ page 157).
.	Volet notification Stocke de l'information sur tous les avertissements et toutes les erreurs des modules du système et tous les événements du contrôleur du système depuis le dernier démarrage du contrôleur du système Pour obtenir des informations complémentaires, reportez-vous à la section Messages () page 356).
\$ _0	Volet paramètres Pour obtenir des informations complémentaires, reportez-vous à la section Volet Paramètres () page 157).
()	Panneau d'information Affiche différentes informations, par exemple, les modules du système attachés et leur numéro de série
M	Panneau de graphiques en ligne (disponible sur le volet accueil) Ouvre une boîte de dialogue indiquant les modules connectés sur la partie supérieure et la valeur physique associée dans la partie inférieure. Le graphique est mis à jour de manière continue et indique le signal des 10 dernières minutes.
∇	Filtre (uniquement disponible sur le volet script et le volet notification) Vous permet de filtrer par module du système, par exemple.

6.6.4 État opérationnel du système

Si le système est connecté à un système de données chromatographiques, l'état opérationnel (par exemple, en veille ou en cours de rinçage) du système apparaît dans le coin supérieur droit de l'écran.

6.6.5 Volet Accueil

Dans le volet accueil, chaque module du système est représenté par une boîte avec une icône spécifique contenant :

- des paramètres fonctionnels, par exemple, la pression de la pompe ;
- une barre colorée en bas de chaque case, représentant la barre à LED. Pour une description des couleurs de la barre à LED, reportezvous à la section Barres à LED (> page 153).



Illustration 84: Boîte de module du système (ici : pompe)

6.6.6 Volet Script

Le volet script offre des scripts automatisés pour plusieurs phases du cycle de vie du produit, par exemple, pour le démarrage et le dépannage.

Vous pouvez sélectionner les scripts pour ouvrir une boîte de dialogue et obtenir de l'information et des instructions détaillées.

6.6.7 Volet Paramètres

Dans ce volet, vous pouvez effectuer les réglages suivants :

- Configurer la Vanquish User Interface :
 - Langue
 - Unités
 - Fuseau horaire
 - Mode de connectivité

- Afficher :
 - le nom du système, la version du microprogramme et les horaires de fonctionnement ;
 - les licences d'utilisation.

6.7 Réglage du dispositif d'affichage

6.7.1 Positions possibles du dispositif d'affichage

Le dispositif d'affichage peut être positionné de quatre manières différentes :

- Position frontale
 - Incliné vers l'arrière



6.7.2 Déplacement du dispositif d'affichage de la position frontale à la position latérale

1. Faites pivoter le dispositif d'affichage de 180 degrés vers la gauche.



Illustration 85: Pivotement du dispositif d'affichage de 180 degrés vers la gauche

2. Inclinez le dispositif d'affichage vers le bas.



Illustration 86: Inclinaison du dispositif d'affichage vers le bas

6.8 Préparer le système à l'utilisation

La présente section contient de l'information sur les étapes supplémentaires requises pour préparer le système à l'utilisation.

- 1. Rincez abondamment le système. Reportez-vous à la section Rinçage du système (▶ page 161).
- 2. Équilibrez le système. Reportez-vous à la section Équilibrage du système (▶ page 163).
- 3. Chargez le carrousel. Reportez-vous à la section Chargement du carrousel (▶ page 165).
- Respectez les recommandations sur les bouchons d'échantillons et les couvercles des plaques à puits. Reportez-vous à la section Recommandations sur les bouchons de flacons et les couvercles des plaques à puits (▶ page 167).

6.8.1 Rinçage du système

AVIS—Les particules et les bulles d'air peuvent nuire au fonctionnement du système

Certains composants du système sont remplis d'isopropanol quand ils sont entreposés.

- Pour éliminer les bulles d'air (et l'isopropanol) avant de redémarrer le système ou après l'installation de composants supplémentaires sur le trajet d'écoulement, exécutez le script : A02 – Démarrer automatiquement.
- Pour éliminer l'isopropanol, utilisez des solvants miscibles à l'isopropanol. S'ils ne le sont pas, utilisez un solvant intermédiaire approprié.

À quel moment exécuter cette opération ?

- Après avoir mis le système en marche
- Après avoir installé de nouveaux composants sur le trajet d'écoulement du système

Liquide requis	Consommation pendant l'exécution du script : A02 – Démarrer automatiquement	Consommation pendant l'analyse	Exigences de liquide
Solvant	Max. 10 mL par réservoir de solvant	En fonction du débit	Lorsque vous utilisez un composant pour la première fois, employez des solvants miscibles à l'isopropanol. S'ils ne le sont pas, utilisez un solvant intermédiaire approprié. Pour tous les autres cas, reportez-vous à la section Informations sur les solvants et les additifs () page 144).
Liquide de rinçage des joints	Max. 10 mL	Environ 250 mL tous les 14 jours	Reportez-vous à la section Remplissage du réservoir de liquide de rinçage des joints () page 134).

Quantités requises de solvant et de liquide de rinçage des joints

Quantités requises de liquide de rinçage de l'aiguille

Consommation pendant le script : A02 – Démarrer automatiquement	Consommatio l'analyse	n pendant	Exigences de liquide
Max. 10 mL de rinçage de l'a injection ¹ :		n par réservoir l'aiguille et par de rinçage de l'aiguille	
	SWP	500 μL	(▶ page 131).
	WWP	400 µL	
	W	250 μL	
¹ : Les valeurs s'appliquent à : boucle d'échantillonnage de 25 µL, injection directe			

Les valeurs s'appliquent a : boucle d'échantillonnage de 25 μL, injection directe (boucle en bypass) configurée dans le système de gestion de données chromatographiques, réglages de rinçage par défaut du système de gestion de données chromatographiques.

Remarque : la consommation dépend de la taille de la boucle d'échantillonnage, la taille de la colonne, la durée du rinçage, le flux du travail et la vitesse de rinçage. Si vous utilisez d'autres réglages, les valeurs pourraient être significativement plus élevées.

Procédez comme suit

- 1. Vérifiez le niveau de liquide de tous les réservoirs, surtout celui des réservoirs de rinçage de l'aiguille. Assurez-vous que les quantités de liquide sont suffisantes pour les points suivants :
 - le script A02 Démarrer automatiquement (reportez-vous au tableau ci-dessus);
 - la première analyse (reportez-vous au tableau ci-dessus et ajoutez les quantités selon le débit de l'application).
- 2. Fermez les portes de tous les modules du système Vanquish Neo, si ce n'est pas déjà fait.
- 3. *Si vous voulez changer de type de solvant* : : pour configurer le type de solvant A/B, exécutez le script : **A01 Définir le type de solvant de la pompe**.
- 4. Cette étape dépend de la situation :
 - Si vous utilisez la colonne installée précédemment : vérifiez et, si nécessaire, ajustez la configuration de la colonne en utilisant les scripts : A03 – Définir le type de colonne de séparation et A04 – Définir les caractéristiques de la colonne de séparation.
 - Si vous avez installé une nouvelle colonne qui n'a pas été utilisée auparavant : configurez la nouvelle colonne à l'aide des scripts :
 A03 – Définir le type de colonne de séparation et A04 – Définir les caractéristiques de la colonne de séparation.
- Exécutez le script avec ou sans diagnostic (test de fuite) : A02 Démarrer automatiquement.
 Le script rince le trajet d'écoulement du système jusqu'au capillaire d'entrée de la colonne, teste la contre-pression du système et effectue un test de fuite, le cas échéant.

6.8.2 Équilibrage du système

À quel moment exécuter cette opération ?

- Après avoir mis le système sous tension et après avoir préparé le système pour l'utilisation initiale
- Après un arrêt de courte durée
- Après avoir changé le débit
- Après avoir changé les températures de consigne

Procédez comme suit

L'équilibrage du système doit comprendre les étapes suivantes :

 Chauffage (ou refroidissement) de tous les appareils du système équipés d'une régulation en température pendant un certain temps :

Étapes d'équilibrage	Durée
Mise à la température initiale (chauffage ou refroidissement) de tous les appareils du système équipés d'une régulation en température. Ce sont, par exemple :	5 heures en fonction des différences de température requises pour l'application
 le débitmètre de pompe ; 	
 la thermostatisation du compartiment à échantillons du passeur d'échantillon ; 	
 le compartiment à colonnes et la colonne. 	

- 2. Pendant le chauffage (ou le refroidissement), effectuez les opérations suivantes selon la situation :
 - Si vous utilisez la colonne installée précédemment : équilibrez la colonne en effectuant le script : B02 – Nettoyer et équilibrer la / les colonne(s).
 - Si vous avez installé une nouvelle colonne qui n'a pas été utilisée auparavant : pour préparer les nouvelles colonnes, exécutez le script : B06 – Conditionnement de la / des colonne(s).
- Surveillez la dérive de la pression de la pompe et vérifiez qu'elle se trouve dans des limites raisonnables pour l'application pendant 5 à 24 heures, selon l'exactitude en matière de temps de rétention nécessaire pour votre application.

6.8.3 Chargement du carrousel

Le carrousel peut accueillir jusqu'à quatre portoirs d'échantillons ou plaques à puits.

Éléments nécessaires

Flacons, portoirs d'échantillons et/ou plaques à puits

Respectez les recommandations de la section Recommandations sur les bouchons de flacons et les couvercles des plaques à puits () page 167).

Préparatifs

- 1. Placez les flacons avec vos échantillons dans le portoir d'échantillon.
- Vérifiez le réglage de la température du compartiment à échantillons et fixez-la tel que requis. Reportez-vous à la section Thermostatage du compartiment du passeur d'échantillon (> page 175).
- 3. Vérifiez la barre à LED du passeur d'échantillon. N'ouvrez pas les portes avant lorsque la barre à LED s'illumine en bleu. Certaines pièces du passeur d'échantillon sont en mouvement.

Procédez comme suit

- Faites pivoter le carrousel afin que le segment souhaité se trouve à l'avant. Vous pouvez faire pivoter le carrousel de la manière suivante :
 - Sélectionnez le bouton ROTATE sur le panneau de commande. Appuyer sur ce bouton fait tourner le carrousel dans le sens inverse des aiguilles d'une montre jusqu'à la position de chargement suivante.
 - Faites pivoter le carrousel manuellement dans le sens désiré.
 - Faites tourner le carrousel pour l'éloigner du ePanel Vanquish Neo dans le système de gestion de données chromatographiques ou dans la case du passeur d'échantillon dans le volet accueil de l'Vanquish User Interface.

 Placez le portoir d'échantillons ou la plaque à puits dans le segment sélectionné. Le portoir d'échantillons ou la plaque à puits doit être bien installé dans le cadre d'alignement et reposer sur les points d'alignement du segment.

AVIS—Orientation correcte des portoirs d'échantillons

Le fait de positionner les portoirs d'échantillons ou les plaques à puits dans le mauvais sens peut endommager le passeur d'échantillon ou donner des résultats erronés.

Positionnez toujours les portoirs d'échantillons et plaques à puits dans le bon sens dans le carrousel, avec la position A1 en haut à gauche.



Illustration 87: Orientation correcte dans le carrousel (ici avec un portoir d'échantillons)

- 3. Chargez les porte-flacons supplémentaires, le cas échéant.
- Faites pivoter le carrousel jusqu'à la position de chargement suivante et chargez-la comme requis. Poursuivez jusqu'à avoir chargé tous les segments requis.
- Fermez les portes avant du passeur d'échantillon. Le passeur d'échantillon démarre un inventaire des portoirs d'échantillons et des récipients d'échantillonnage présents dans le compartiment à échantillons.

Identification automatique du type de portoir

Si vous utilisez des portoirs d'échantillons ou des plaques à puits avec code-barres Vanquish, le type et l'orientation des portoirs sont identifiés automatiquement.

Configuration manuelle du type de portoir

Si vous utilisez des portoirs d'échantillon ou des plaques à puits sans code-barres Vanquish, vous devez configurer manuellement le type de portoir de chaque segment du ePanel de Vanquish Neo dans le système de gestion de données chromatographiques.

Consultez également

- Carrousel () page 50)
- Identification du type de portoir (> page 51)

6.8.4 Recommandations sur les bouchons de flacons et les couvercles des plaques à puits

Utilisez seulement des bouchons de flacons et des couvercles de plaques à puits selon les recommandations ci-dessous.

Bouchons de flacons et septum

Utilisez seulement du septum sans talc à base de silicone non rempli/ PTFE. Les particules de talc ou de remplissage des joints pourraient causer des problèmes de contre-pression ou de blocages dans le système.

Exemple de bouchons répondant aux exigences : bouchons de flacons à visser et septum collés Thermo Scientific en PTFE rouge/silicone blanc

Tapis de couverture pour plaques à puits

Utilisez des tapis de couverture non remplis en silicone.

Exemple de tapis répondant aux exigences : tapis en silicone Thermo Scientific WebSeal[™] microplaque non revêtue à 96 puits en plastique WebSeal

Ruban d'étanchéité pour plaque à puits

Utilisez du ruban d'étanchéité et de fermeture *sans* adhésif sur la surface des cavités. La présence d'adhésif sur cette surface pourrait contaminer l'aiguille d'injection.

Exemple de ruban répondant aux exigences : Thermo Scientific WebSeal microplaque non revêtue à 96 puits en plastique avec ruban d'étanchéité WebSeal avec 3 couches sans adhésif à la surface des cavités.

N'utilisez *pas* des joints de plaque à puits susceptibles de faire en sorte que le matériau du joint soit extrait par l'aiguille du passeur d'échantillon. Par exemple, des parties d'un ruban en aluminium pourraient être extraites d'un ruban d'étanchéité pour plaques à puits en aluminium et se déposer dans le siège de l'aiguille.

6.9 Optimisation des performances

La présente section fournit des informations sur les performances optimales et vous indique comment les améliorer encore davantage.

6.9.1 Interdépendance entre la pression de la pompe et le débit

La pression maximale que la pompe peut livrer dépend des paramètres suivants :

- Débit sélectionné
- Viscosité du solvant utilisé dans les voies (voie A ou voie B)

Pour les solvants pré-étalonnés, le graphique montre comment la pression de pompe actuelle dépend du débit et de la voie sélectionnés.



Illustration 88: Dépendance de la pression de la pompe et du débit pour la voie A et la voie B

Pour les autres solvants, respectez les consignes suivantes :

- Pour les débits de gradient jusqu'à 10 μL/min, utilisez un solvant aqueux (faible) dans la voie A et un solvant organique (fort) dans la voie B.
- Si le débit dépasse 10 μL/min et la contre-pression attendue du système est supérieure à 140 MPa (1 400 bar), utilisez le solvant présentant la viscosité la plus élevée dans la voie A.

Pendant l'utilisation, la pompe surveille continuellement s'il est possible d'offrir le débit sélectionné à la pression résultante. Si la contre-pression du système devient si élevée que la pompe ne peut offrir le débit sélectionné, voici qui se passe :

- L'injection ou la séquence arrête.
- Un message apparaît indiquant, par exemple, que la pompe ne peut offrir le débit demandé à la pression actuelle (**code 4531**) ou que la pression dépasse la limite absolue (**code 4503**).

Lorsque vous réduisez le débit, la contre-pression effective du système diminue. Dans le même temps, la quantité de pression disponible à partir de la pompe augmente. Si la pompe ne parvient toujours pas à fournir le débit sélectionné, réduisez davantage la contre-pression du système. Par exemple, en réduisant encore le débit, en augmentant la température de la colonne, ou en augmentant le diamètre intérieur des capillaires fluidiques.

6.9.2 Ajout de solvants personnalisés au tableau d'étalonnage

La pompe peut offrir des débits précis seulement si le solvant que vous souhaitez utiliser est étalonné. Les solvants étalonnés peuvent être sélectionnés dans le script **A01 – Définir le type de solvant de la pompe** ou dans le tableau de calibration des solvants. Les types et les mélanges de solvants les plus courants ont été étalonnés à l'usine :

- Eau, 100 %
- Acétonitrile, 100 %
- Mélange d'acétonitrile et d'eau, 80:20 (v:v)
- Mélange de méthanol et d'eau, 90:10 (v:v)

Si vous souhaitez utiliser d'autres solvants, étalonnez-les une fois pour les ajouter à la configuration. L'étalonnage est recommandé pour les types de solvants s'écartant de > 5 % (volume par volume) d'un des solvants mentionnés ci-dessus.

Exemples

Exemple 1

Solvant		Action
A	98:2 (v:v), eau:acétonitrile plus 0,1 % d'acide formique	Utilisez les données d'étalonnage de 100 % d'eau.
В	80:20 (v:v), acétonitrile:eau plus 0,1 % d'acide formique	Utilisez les données d'étalonnage de 80:20 % acétonitrile.

Exemple 2

Solvant		vant	Action
	A	95:5 (v:v), eau:diméthylsulfoxyde plus 0,1 % d'acide formique	Utilisez les données d'étalonnage de 100 % d'eau.
-	В	75:20:5 (v:v:v), acétonitrile:eau:diméthylsulfoxyde plus 0,1 % d'acide formique	Utilisez les données d'étalonnage de 80:20 % acétonitrile.

Exemple 3

Solvant		Action
A	100 % d'eau plus 0,1 % d'acide formique	Utilisez les données d'étalonnage de 100 % d'eau.
В	90:10 (v:v), acétonitrile:eau plus 0,1 % d'acide formique	L'étalonnage est requis. Reportez- vous aux explications plus bas.

Notes générales

- L'étalonnage de solvant est effectué par voie.
- Si vous souhaitez utiliser un solvant dans les deux voies, vous devez effectuer l'étalonnage deux fois pour étalonner séparément le solvant de chaque voie.
- L'étalonnage réel est automatisé et s'exécute sans surveillance. Avant de lancer le script de l'étalonnage réel, vous devez régler et configurer la pompe en vue de l'étalonnage et fournir des entrées sur l'utilisateur. Une fois l'étalonnage terminé, vous devez retourner la pompe à la configuration capillaire d'origine et raccorder tous les capillaires et le système selon les besoins de l'application.
- Les scripts d'étalonnage vous guident à travers toutes les étapes du processus.
- Le temps requis pour achever l'étalonnage réel est environ 100 minutes. De plus, prévoyez 30 minutes avant pour préparer la configuration de l'étalonnage et 30 minutes après pour restaurer la configuration opérationnelle.

Pièces et éléments nécessaires

- Solvant convenant à la purge, afin d'évacuer toute substance nocive
- Solvant à étalonner
- Bouchon de raccordement Viper (3 bouchons)

Procédez comme suit

- 1. Pour volet script, paramétrez le filtre sur Étalonnage du solvant.
- 2. Exécutez et suivez les scripts d'étalonnage de solvant l'un après l'autre.
- 3. Après avoir rétabli la configuration d'utilisation, vérifiez si la pompe fuit. Exécutez le script : **E04 Test de fuite détaillé (pompe)**.

Résultats d'étalonnage

- Étalonnage réussi : le solvant est ajouté au tableau d'étalonnage de solvants et vous pouvez maintenant le sélectionner.
- Une erreur est survenue pendant l'étalonnage : à partir des scripts d'étalonnage, complétez l'étalonnage ou rétablissez la configuration d'utilisation.

6.9.3 Changement de type de solvant

- 1. Si vous souhaitez changer de type de solvant :
 - a) Pour configurer le type de solvant A/B, exécutez le script : **A01 – Définir le type de solvant de la pompe**.
 - b) Pour purger le système, exécutez le script : B01 Modifier les liquides / solvants (avec les paramètres suivants : Solvant A/B et Rafraîchir Uniquement = désactivé).

6.9.4 Changement du liquide de rinçage des joints

Si vous souhaitez changer de liquide de rinçage des joints :

Exécutez le script : **B01 – Modifier les liquides / solvants** (avec les paramètres suivants : **Liquide de rinçage des joints arrière** et **Rafraîchir Uniquement = désactivé**).

6.9.5 Changement du liquide de rinçage de l'aiguille

Si vous souhaitez changer de liquide de rinçage de l'aiguille :

Exécutez le script : **B01 – Modifier les liquides / solvants** avec les paramètres de rinçage de l'aiguille respectifs et **Rafraîchir Uniquement = désactivé**.

6.9.6 Réglage des décalages du capteur de débit de la pompe

Le script de réglage du décalage du capteur de débit mesure le signal du capteur sans débit et enregistre les résultats pour qu'il soit possible de corriger le signal de flux en conséquence. Ainsi, le changement de décalage qui s'est produit depuis l'étalonnage à l'usine sera supprimé.

Réglage non requis

Il n'est pas nécessaire de régler les décalages du capteur de débit quand les temps de rétention sont stables ou à l'intérieur des limites attendues. Ne réglez pas les décalages du capteur de débit dans les situations suivantes :

- Pendant une étude scientifique ou une série de mesures.
- Pendant une série de mesures qui doit donner des résultats comparables.
- Très fréquemment, par exemple tous les jours.

Réglage potentiellement requis

Afin d'obtenir des résultats de mesures précis et comparables, il peut être nécessaire de régler les décalages du capteur de débit dans les situations suivantes :

- Quand les temps de rétention sont significativement différents des limites attendues (décalés). Cette situation peut entraîner des pics ratés.
- Quand les temps de rétention varient considérablement comparativement aux mesures précédentes.

Il est estimé que le réglage des décalages du capteur de débit sera requis une fois tous les mois ou tous les trois mois, surtout en effectuant des analyses < 2 μ L/min.

Procédez comme suit

Respectez les consignes suivantes avant de commencer :

- Le système doit être allumé depuis au moins 5 heures. Seulement à ce moment, un réglage sera possible.
- *Recommandation* : effectuez le réglage en utilisant les solvants de votre application.
- Exécutez le script de réglage suivant : C01 Régler les décalages du capteur de débit de la pompe.
- Quand le réglage est terminé, prévoyez un temps d'équilibrage pour obtenir une stabilité optimale du temps de rétention, par exemple, un cycle à vide.

6.9.7 Rinçage de l'aiguille

Le cycle de rinçage de l'aiguille rince l'aiguille pour en éliminer tout échantillon résiduel. Ce cycle peut être compris dans la séquence de prélèvement des échantillons ou réalisé manuellement. Cette section décrit comment lancer un cycle de rinçage manuel de l'aiguille pour optimiser la performance du système et de la séquence du flux de travail.

À quel moment exécuter cette opération ?

- Remplacement du liquide de rinçage de l'aiguille
- Remplacement des conduites de rinçage
- Maintenance préventive et de routine pour le dispositif de rinçage de l'aiguille

Préparatifs

Respectez les consignes relatives au rinçage de l'aiguille, reportez vous à la section Remplissage des réservoirs de rinçage de l'aiguille () page 131).

Procédez comme suit

Pour commencer le rinçage extérieur de l'aiguille, suivez les étapes suivantes :

Appuyez sur le bouton **WASH** sur le panneau de commande.

Cela lance un cycle de rinçage extérieur de l'aiguille au cours duquel la surface externe de l'aiguille est rincée dans l'orifice de rinçage.

Pour commencer le rinçage manuel intérieur et extérieur de l'aiguille, suivez les étapes suivantes :

Sur le VSC, exécutez le script : **B03 – Nettoyer le système** Ainsi, le système est initialisé pour effectuer une procédure de rinçage de l'extérieur et de l'intérieur de l'aiguille, y compris la boucle d'échantillonnage.

Une fois le cycle de rinçage terminé, l'aiguille retourne dans son siège.

Consultez également

- Dispositif de rinçage de l'aiguille (vue d'ensemble) (> page 58)
- Entretien du dispositif de rinçage de l'aiguille () page 273)

6.9.8 Thermostatage du compartiment du passeur d'échantillon

Dans le passeur d'échantillon, vous pouvez effectuer un thermostatage du compartiment.

Procédez comme suit

- Activez la régulation de la température dans le système de gestion de données chromatographiques ou dans la case du passeur d'échantillon dans le volet accueil de l'Vanquish User Interface.
- 2. Réglez la température cible. Vérifiez que la température de thermostatage convient à tous les échantillons.

NOTE Si vous arrêtez le passeur d'échantillon, il conserve la température configurée. Vérifiez la température quand vous le rallumez et adaptezla, le cas échéant.

6.9.9 Thermostatage du compartiment à colonnes

Vous pouvez effectuer un thermostatage du compartiment à colonnes.

Procédez comme suit

- 1. Activez la régulation de la température dans l'Vanquish User Interface ou en appuyant sur le bouton **TEMP CTRL**.
- Réglez la température cible dans le système de gestion de données chromatographiques ou dans la case du compartiment à colonnes dans le volet accueil de l'Vanquish User Interface. Vérifiez que la température de thermostatage convient aux pièces suivantes :
 - Vannes
 - Colonnes

Notes sur le réglage de la température

Le compartiment à colonnes n'est pas muni d'un système de refroidissement actif. Veuillez noter les points suivants :

- Vous ne pouvez pas diminuer la température du compartiment à colonnes en dessous de la température venant de la température ambiante et de l'auto-échauffement du module.
- La température réglée doit être au moins 5°C au-dessus de la température ambiante.

Ready State			
	La propriété Ready du système de gestion de données chromatographiques indique si le compartiment à colonnes est prêt à fonctionner. Quand l'état du compartiment à colonnes est NotReady , vous ne pouvez pas lancer l'analyse. Les paramètres Ready Temp Delta et Equilibration Time influencent la durée nécessaire au compartiment à colonnes pour passer à l'état Ready .		
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'équilibrag	age		
	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. 		

6.10 Arrêt du système

Si une période d'inactivité du système est prévue, suivez les instructions de cette section.

6.10.1 Interruption du fonctionnement pour une courte durée

Pour interrompre le fonctionnement du système pendant une courte durée (arrêt de courte durée), par exemple la nuit, suivez les instructions suivantes :

Préparatifs

Si la configuration de votre système comprend un compartiment à colonnes, réglez la température maximale à 40°C.

Procédez comme suit

1. Exécutez le script et suivez les instructions : **B05 – Arrêt (de courte durée)**.

Redémarrage du système après un arrêt à court terme

- 1. Avant de reprendre l'utilisation, suivez les étapes suivantes :
 - a) Laissez le débit de la pompe et la température s'équilibrer pour retourner aux conditions d'utilisation.
 - b) Purgez le passeur d'échantillon avec du liquide de rinçage de l'aiguille faible. Exécutez le script : B01 – Modifier les liquides / solvants avec les paramètres Liquide de rinçage de l'aiguille faible pour rinçage interne et Rafraîchir Uniquement = activé.
- 2. Exécutez le script : B02 Nettoyer et équilibrer la / les colonne(s).

6.10.2 Interruption du fonctionnement pour une longue période

Pour interrompre le fonctionnement pour une période prolongée et garder le système dans le laboratoire, suivez les instructions ci-dessous.

 Exécutez le script : B05 – Arrêt (de longue durée) avec le paramètre suivant : Module à retirer = aucun.

NOTE Il n'est généralement pas nécessaire de retirer les conduites de solvant des guides de tubulure (même pour remplacer le module à glissière). Toutefois, lorsque vous retirez les conduites de solvant des guides de tubulure dans la pompe et dans tous les modules placés audessus de la pompe dans le système empilé, veillez à ne pas tirer sur les autres tubes présents dans les guides.

- 2. Videz les réservoirs de solvant.
- 3. Arrêtez le système au moyen du bouton d'alimentation électrique (bouton du bas sur le socle du système).
- 4. Arrêtez le contrôleur du système au moyen de son bouton d'alimentation électrique (bouton du haut sur le socle du système).

NOTE Lorsque le passeur d'échantillon est hors tension, sa porte avant gauche s'ouvre automatiquement afin que le compartiment à échantillons soit correctement ventilé et elle ne peut pas être fermée.

Redémarrage du système après un arrêt à long terme

- Remplissez les réservoirs de solvant. Reportez-vous à la section Remplissage des réservoirs de solvant (▶ page 129).
- 2. Mettez le système sous tension. Reportez-vous à la section Première mise en marche du système (▶ page 136).
- 3. Préparez le système à l'utilisation. Reportez-vous à la section Préparer le système à l'utilisation (▶ page 161).

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.

Les procédures de maintenance ne nécessitent pas de retirer les portes. 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 des portes (> page 333).
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 26).



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—Aiguille d'injection à pointe acérée

L'aiguille d'injection est dotée d'une pointe très acérée qui peut causer des blessures cutanées.

Pour éviter les blessures, ne touchez jamais la pointe de l'aiguille.



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 du système 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 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. Reportez-vous à la section Vidange des conduites de solvant (> page 237).
- 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.

AVIS

Déplacer le système empilé Vanquish, même sur une courte distance sur la paillasse de laboratoire, peut endommager le socle du système et altérer son fonctionnement.

Avant de déplacer le système empilé, déverrouillez le socle du système à l'aide des outils de déverrouillage.

NOTE Lorsque le passeur d'échantillon est hors tension, sa porte avant gauche s'ouvre automatiquement afin que le compartiment à échantillons soit correctement ventilé, et elle ne peut pas être fermée tant que le passeur d'échantillon est hors tension.

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.

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.
- Après les travaux de maintenance et d'entretien sur les raccordements fluidiques du système, effectuez les étapes suivantes : premièrement, testez le système pour vérifier la contrepression (script : D01 – Tester la contre-pression du système), puis vérifiez que le système ne présente aucune fuite (script : D02 – Tester le serrage du système (ce qui devrait être testé = le système)).

AVIS

Les raccordements fluidiques et raccords de capillaires sont extrêmement sensibles à la contamination. La poussière et les débris peuvent contaminer ces raccordements.

Installez toujours des capuchons sur les capillaires et des bouchons sur les raccordements fluidiques ouverts pour éviter toute contamination.

Réparations en atelier

• 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.

Consultez également

Consommables et pièces de rechange (page 417)

7.4 Maintenance de routine et maintenance préventive

Seul un système 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 les tableaux de maintenance suivants. Les fréquences indiquées dans les tableaux 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 le système.

7.4.1.1 Maintenance hebdomadaire

Composant du produit	Opération de maintenance à réaliser
Système/tous les modules	 Raccordements fluidiques : Inspectez les raccordements fluidiques afin d'y détecter d'éventuels signes de :
	 Blocage : Exécutez le script et suivez les instructions : D01 – Tester la contre-pression du système.
	 Dépôts de sel
	 Bulles d'air Agitez les solvants et purgez la pompe et le passeur d'échantillon, le cas échéant.
	 Solvants : Si vous utilisez l'un des solvants suivants, procédez comme suit :
	 Eau pure ou un tampon aqueux : Pour remplacer le réservoir de solvant, exécutez le script : B01 – Modifier les liquides / solvants.
	 Tampons ou solutions salines : rincez soigneusement le système après utilisation avec un solvant adéquat ne contenant ni tampon ni sel. Reportez-vous à la section Purge (vue d'ensemble) (▶ page 65).
	 Solvants corrosifs ou méthanol avec < 10 % d'eau : rincez le système avec de l'eau ou d'autres solvants appropriés non corrosifs. Reportez-vous à la section Purge (vue d'ensemble) () page 65).

Composant du produit	Opération de maintenance à réaliser
Dispositifs de rinçage	 Vérifiez le niveau de liquide. Vérifiez si les réservoirs de rinçage contiennent des particules, de la poussière ou des algues. Si c'est le cas, remplacez les réservoirs de rinçage par des réservoirs propres. Exécutez le script : B01 – Modifier les liquides / solvants (Rafraîchir Uniquement = activé). Assurez-vous que les réservoirs de rinçage ne contiennent ni détergents ni substances chimiques résiduels. Nous recommandons de rincer les réservoirs de rinçage avec une petite quantité du liquide de rinçage qui sera utilisé dans le système.
Passeur d'échantillon	 Inspectez les échantillons et plaques à puits dans le passeur d'échantillon afin d'y détecter d'éventuels fissures ou défauts. Nettoyez les déversements dans le passeur d'échantillon, le cas échéant.

7.4.1.2 Maintenance régulière

Composant du produit	Opération de maintenance à réaliser
Système/tous les modules	 Remplacez les réservoirs de solvant toutes les semaines ou toutes les deux semaines. Pour les remplacer, exécutez le script : B01 – Modifier les liquides / solvants. Avant de remplir les réservoirs, rincez-les soigneusement.
	 Vérifiez si le système fuit : D02 – Tester le serrage du système (ce qui devrait être testé = le système).
	 Remplacez le filtre de conduite de solvant à des intervalles réguliers. Ceci est particulièrement important si vous utilisez des solvants aqueux. Les solvants aqueux peuvent contaminer les filtres en raison de la formation et du dépôt d'algues et de microorganismes sur les frittés des filtres.
	 Vérifiez que les raccordements fluidiques sont en bon état. Recherchez des fissures, des entailles, des coupures et des obstructions.
	 Assurez-vous que le dispositif d'évacuation n'est pas obstrué. Reportez-vous à la section Vérification du dispositif d'évacuation (> page 128).
	 Videz les bacs d'évacuation.
	 Nettoyez le système. Reportez-vous à la section Nettoyage ou décontamination d'un module du système ou du système (> page 191).
	 Inspectez les étiquettes d'avertissement sur le système et les modules. Si une étiquette d'avertissement est manquante ou illisible, contactez Thermo Fisher Scientific pour obtenir une étiquette de remplacement.

Composant du produit	Opération de maintenance à réaliser
Dispositifs de rinçage	 Dispositifs de rinçage de l'aiguille : Remplacez régulièrement les réservoirs de rinçage de l'aiguille, environ toutes les semaines ou toutes les deux semaines. Pour remplacer ces réservoirs, exécutez le script : B01 – Modifier les liquides / solvants.
	 Système de rinçage des joints : Remplacez le réservoir de liquide de rinçage des joints. Rincez le réservoir soigneusement avant tout remplissage. Utilisez un solvant ultrapur pour le rinçage. Inspectez les conduites de rinçage des joints. Si une conduite est obstruée (pincée ou pliée), remplacez-la. Inspectez le tube péristaltique de la pompe de rinçage des joints. S'il est déplacé ou durci, replacez ou remplacez le tube. Reportez-vous à la section Remplacement du tube péristaltique dans la pompe de rinçage des joints (> page 293). Remplacez les tubes tous les 6 mois à titre de maintenance préventive. Vérifiez s'il y a des fuites dans le système de rinçage des joints. Reportez-vous à la section Test d'étanchéité du système de rinçage des joints (> page 282).
Pompe	 Nettoyez les clapets de non-retour de la tête de pompe une fois par mois. Reportez-vous à la section Nettoyage des clapets de non-retour (tête de pompe) (> page 229).
	 Rincez la pompe avec des solvants fraîchement préparés pour éviter toute contamination. Reportez-vous à la section Rinçage de la pompe pour éviter toute contamination (> page 189).
Passeur d'échantillon	 Remplacez les tubes de la pompe d'évacuation tous les 6 mois. Reportez-vous à la section Remplacement du tube de la pompe d'évacuation () page 270). (Chauffé) Flux de tâches Retenir-et-Éluer uniquement : déterminez la contre-pression du trajet d'écoulement de la colonne de piégeage. Exécutez le script : D01 – Tester la contre- pression du système avec le paramètre : Tester uniquement la résistance de la colonne de piégeage. Prenez note de la valeur et comparez-la aux tests de contre-pression ultérieurs pour évaluer le vieillissement de la colonne.
Dispositif d'affichage	Resserrez les vis qui fixent le dispositif d'affichage au système.

Consultez également

Utilisation de solvants et d'additifs (> page 145)

7.4.1.3 Maintenance annuelle

Composant du produit	Opération de maintenance à réaliser
Système/tous les modules	Demandez au personnel d'entretien de Thermo Fisher Scientific d'effectuer une maintenance préventive une fois par an.
Système de rinçage des joints	 Remplacez les joints de rinçage des joints. Reportez-vous à la section Remplacement du joint de rinçage des joints (> page 216).
	 Remplacez les conduites de rinçage des joints. Reportez-vous à la section Entretien du système de rinçage des joints (> page 282).
Pompe	Inspectez les pistons à la recherche de dépôts de particules ou d'éventuels signes de détérioration, stries ou rayures. Reportez- vous à la section Nettoyage du piston () page 203).

7.4.2 Rinçage de la pompe pour éviter toute contamination

Matériel nécessaire

Selon les besoins :

- Solvant fraîchement préparé
- Isopropanol
- Éthanol

Respectez les consignes suivantes

Les consignes suivantes s'appliquent à l'utilisation classique du système Vanquish recommandée dans le présent manuel, avec notamment l'emploi de solvants fraîchement préparés de haute qualité (au moins de qualité UHPLC ou LC/MS) et fréquemment remplacés :

- Il suffit généralement de rincer la pompe avec des solvants fraîchement préparés.
- Si vous utilisez de l'eau, de l'acétonitrile ou du méthanol pour la chromatographie, il convient de rincer la pompe une fois par semaine.
- Les phases mobiles contenant des sels, des tampons et/ou d'autres additifs sont par nature sujettes à la contamination organique et/ou à la biodégradation. Quand vous utilisez des tampons non volatils, respectez les consignes suivantes :
 - filtrez les phases mobiles avant utilisation ;
 - rincez le système avec, par exemple, de l'isopropanol > 60 % ou de l'éthanol toutes les semaines pour pouvoir continuer à l'utiliser sans problème.

• Adaptez cet intervalle en fonction des solvants utilisés.

NOTE En cas de contamination persistante, par exemple si des pics fantômes reproductibles apparaissent sur le chromatogramme vierge sans injection d'échantillon, envisagez une procédure de nettoyage plus rigoureuse. Reportez-vous à la section Rinçage de la pompe en cas de contamination persistante uniquement (> page 190).

Consultez également

Utilisation de solvants et d'additifs (> page 145)

7.4.3 Rinçage de la pompe en cas de contamination persistante uniquement

À quel moment exécuter cette opération ?

NOTE Si le système Vanquish Neo est utilisé de manière classique, comme recommandé dans le présent manuel, il suffit généralement de rincer la pompe régulièrement. Reportez-vous à la section Rinçage de la pompe pour éviter toute contamination () page 189).

Par exemple, si des pics fantômes apparaissent de façon répétée sur le chromatogramme vierge sans qu'il y ait injection d'échantillon, envisagez d'utiliser une procédure de nettoyage plus rigoureuse. Suivez les instructions ci-dessous.

Pièces et éléments complémentaires nécessaires

- Capillaire de contre-pression adapté, générant environ 20 à 90 MPa à un débit de 50 μl/min dans l'eau (par exemple, capillaire nanoViper, verre de silice, diamètre intérieur : 20 μm, longueur : 350 mm)
- 6 moles (6 M) d'acide nitrique (qualité HPLC)
- Bac d'évacuation (pour collecter l'acide nitrique)
- Eau fraîche (qualité HPLC)
- Acétonitrile fraîchement préparé (qualité HPLC)
- Nouveaux réservoirs de solvant avec solvant fraîchement préparé
- Nouveaux frittés pour filtres de conduite de solvant

Procédez comme suit

 Installez le capillaire de contre-pression sur l'orifice de refoulement de la pompe. Dirigez l'extrémité libre du capillaire vers le bac d'évacuation.

- 2. Remplacez les frittés dans les filtres de conduite de solvant par de nouveaux frittés.
- Remplacez le bac d'évacuation installé par le nouveau bac d'évacuation pour éviter toute réaction chimique entre l'acide nitrique et le solvant dans le bac d'évacuation.
- 4. Rincez la pompe avec 6 M d'acide nitrique pendant 2 heures, en appliquant 50 % de B à un débit de 50 μ L/min.
- 5. Rincez la pompe avec de l'eau fraîche de qualité HPLC jusqu'à ce que le pH soit neutre.
- 6. Retirez le bac d'évacuation contenant l'acide nitrique et réinstallez le bac d'évacuation précédemment installé.
- 7. Rincez la pompe avec de l'acétonitrile fraîchement préparé pendant 2 heures, en appliquant 50 % de B à un débit de 50 μ L/min.
- 8. Raccordez les nouveaux réservoirs de solvant contenant le solvant fraîchement préparé.
- 9. Désinstallez le capillaire de contre-pression et raccordez le système selon les besoins de votre application.
- 10. Purgez la pompe.
- 11. Équilibrez le système.

Consultez également

Purge (vue d'ensemble) (▶ page 65)

7.4.4 Nettoyage ou décontamination d'un module du système ou du système

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

La décontamination est requise, par exemple, en cas de fuite ou de déversement accidentel, ou avant l'entretien ou le transport de la pompe. Utilisez un détergent ou désinfectant approprié afin de vous assurer que le traitement permette une manipulation sans danger du système et de ses modules. 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 avec de 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. Si vous souhaitez nettoyer ou décontaminer seulement le passeur d'échantillon ou le système :
 - a) Déplacez l'unité d'aiguille à la position d'entretien.
 - b) Resserrez le vis de verrouillage d'expédition (reportez-vous à la section Verrouillage de l'unité d'aiguille (▶ page 251)).
 - c) Retirez tous les récipients d'échantillon et les portoirs d'échantillons du compartiment à échantillons.
- 2. Procédez comme suit :
 - Si vous souhaitez nettoyer ou décontaminer un module du système en particulier : Mettez le module du système hors tension et débranchez le cordon d'alimentation de la source d'alimentation.
 - Si vous souhaitez nettoyer ou décontaminer le system : Mettez tous les modules du système et le système hors tension et débranchez les cordons d'alimentations de la source d'alimentation.

Procédez comme suit

- 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.

Étapes à suivre après le nettoyage ou la décontamination

- 1. Si vous avez nettoyé ou décontaminé seulement le passeur d'échantillon ou le système :
 - a) Desserrez la vis de verrouillage d'expédition.
 - b) Placez les récipients d'échantillonnage et les portoirs d'échantillons dans le compartiment à échantillons.
- 2. Branchez le(s) cordon(s) d'alimentation sur la source d'alimentation.

- Mettez le(s) module(s) du système sous tension.
 Après le démarrage du passeur d'échantillon, l'unité d'aiguille se déplace automatiquement à la position Inject.
- 4. Préparez le système pour utilisation (reportez-vous à la section Préparer le système à l'utilisation (▶ page 161)).

7.5 Maintenance de la tête de pompe

La présente section décrit comment entretenir ou remplacer la tête de pompe ou ses pièces. Suivez les instructions de la procédure de maintenance que vous souhaitez effectuer.

- Pour remplacer une tête de pompe, reportez-vous à la section Remplacement de la tête de pompe (▶ page 197).
- Pour remplacer un piston, reportez-vous à la section Remplacement des pistons (> page 201).
- Pour nettoyer un piston, reportez-vous à la section Nettoyage du piston (▶ page 203).
- Pour remplacer une bague de support, reportez-vous à la section Remplacement de la bague de support (▶ page 204).
- Pour remplacer un joint de piston, reportez-vous à la section Remplacement du joint du piston (▶ page 209).
- Pour remplacer un joints de rinçage des joints, reportez-vous à la section Remplacement du joint de rinçage des joints (> page 216).
- Pour remplacer un joint de la tête de pompe, reportez-vous à la section Remplacement des joints de la tête de pompe (> page 220).
- Pour nettoyer ou remplacer un clapet de non-retour, reportez-vous à la section Entretien des clapets de non-retour de la tête de pompe (> page 226).

Consultez également

Vue d'ensemble des éléments de la tête de pompe (> page 196)

7.5.1 Vue d'ensemble des éléments de la tête de pompe



L'image montre les principaux éléments de la tête de pompe.

Illustration 89: Vue d'ensemble des éléments de la tête de pompe

N°	Description
1	Tête de pompe
2	Clapet de non-retour d'aspiration
3	Joints de piston
4	Bagues de support
5	Joints de tête de pompe, pour les vis de la plaque de rinçage des joints (joints toriques 6,5x1)
6	Joints de tête de pompe, pour les vis de la plaque de rinçage des joints (joints toriques 11x1,5)
7	Plaques de rinçage des joints
8	Vis, plaques de rinçage des joints
9	Joints de rinçage des joints
10	Pistons
11	Raccords pour tube de rinçage des joints (à visser)
12	Clapet de non-retour de refoulement
13	Tube en U

7.5.2 Remplacement de la tête de pompe

À quel moment exécuter cette opération ?

Détérioration de tête de pompe

Procédez comme suit

- Retirez la tête de pompe. Reportez-vous à la section Retrait de la tête de pompe (▶ page 197).
- 2. Installez la tête de pompe. Reportez-vous à la section Installation de la tête de pompe (▶ page 198).

7.5.2.1 Retrait de la tête de pompe

À quel moment exécuter cette opération ?

- Pour remplacer une tête de pompe
- Pendant la maintenance des éléments de la tête de pompe

Outils et éléments complémentaires nécessaires

- Clé hexagonale, taille 6
- Solvant convenant pour la purge, afin d'évacuer toute substance nocive

Préparatifs

1. Afin d'évacuer toutes les substances nocives, purgez la pompe à l'aide d'un solvant approprié.

Éléments touchés sur la tête de pompe



Illustration 90: Tête de pompe

N°	Description
1	Câble du capteur de pression
2	Conduite de solvant sur le clapet de non-retour d'aspiration

N°	Description
3	Vis de la tête de pompe
4-6	Conduites de rinçage des joints
7	Capillaire raccordant la tête de pompe et le débitmètre

Procédez comme suit

- Pour la tête de pompe sur laquelle vous voulez exécuter la procédure, déconnectez les pistons depuis le panneau de commande (boutons SELECT et DOCK) ou le ePanel dans le système de gestion de données de chromatographie.
- Attendez que les pistons soient déconnectés avant de procéder. La déconnexion est terminée une fois que la pompe s'est arrêtée et que la LED située à côté du bouton FLOW clignote en vert pour la tête de pompe sélectionnée.
- 3. Débranchez le câble du capteur de pression du connecteur **Pressure**.
- 4. Débranchez la conduite de solvant du clapet de non-retour d'aspiration.
- 5. Retirez le capillaire qui raccorde la tête de pompe du débitmètre.
- 6. Retirez les conduites de rinçage des joints des raccords de tube sur la tête de pompe. Ne retirez pas les raccords de tube.
- 7. Desserrez la vis de tête de pompe avec la clé hexagonale.
- 8. Maintenez la tête de pompe d'une main et retirez la vis de la tête de pompe.
- Retirez la tête de pompe en tirant sur la tête d'environ 10 mm (0,5 pouce), en écartant la pompe du panneau avant et en l'inclinant vers le bas.

Les pistons sont généralement retirés de la pompe en même temps que la tête de pompe. Si un piston reste dans la pompe, inclinez-la et retirez le piston avec vos doigts.

Consultez également

- Installation de la tête de pompe (> page 198)
- Purge (vue d'ensemble) (> page 65)

7.5.2.2 Installation de la tête de pompe

À quel moment exécuter cette opération ?

- Pour installer une tête de pompe de remplacement
- Pendant la maintenance des éléments de la tête de pompe

Éléments nécessaires

Tête de pompe

Outils et éléments complémentaires nécessaires

- Clé hexagonale, taille 6
- Solvant convenant à votre application

Éléments touchés sur la tête de pompe



Illustration 91: Tête de pompe

N°	Description
1	Câble du capteur de pression
2	Conduite de solvant sur le clapet de non-retour d'aspiration
3	Vis de la tête de pompe
4-6	Conduites de rinçage des joints
7	Capillaire raccordant la tête de pompe et le débitmètre

Respectez les consignes suivantes

Situation	Description
Réinstallation de la tête de pompe	N'intervertissez pas les têtes de pompe. Prenez note de l'étiquetage sur le câble du capteur de pression (A pour la tête de pompe gauche, B pour la tête de pompe droite).
	Si vous devez réinstaller un piston, reportez-vous à la section Installation des pistons () page 202).
Installation d'une nouvelle tête de pompe (tête de pompe de remplacement)	Lors d'une nouvelle installation, les nouvelles têtes de pompe peuvent être installées dans l'une ou l'autre des positions. Elles ne portent pas d'étiquette sur le câble du capteur de pression. Envisagez d'y apposer des étiquettes indiquant leur position d'installation en prévision de futures opérations de maintenance.

Procédez comme suit

- 1. Vérifiez que les pistons sont insérés aussi loin que possible dans la tête de pompe.
- 2. Insérez la tête de pompe dans la pompe et serrez la vis de tête de pompe à l'aide de la clé hexagonale.
- 3. Raccordez les conduites de rinçage des joints aux raccords de tube sur la tête de pompe.
- 4. Branchez le capillaire qui raccorde la tête de pompe au débitmètre.
- 5. Branchez la conduite de solvant sur le clapet de non-retour d'aspiration.
- 6. Branchez le câble du capteur de pression sur le connecteur **Pressure**.
- Connectez les pistons depuis le panneau de commande (boutons SELECT et DOCK) ou le ePanel dans le système de gestion de données de chromatographie.
- Attendez que les pistons soient connectés avant de procéder. La connexion est terminée une fois que la pompe s'est arrêtée et que la LED située à côté du bouton FLOW arrête de clignoter en vert pour la tête de pompe sélectionnée.
- 9. Mettez la pompe hors tension au moyen de son interrupteur d'alimentation principal, puis rallumez-la.
- 10. Purgez la pompe à l'aide du solvant convenant à votre application.
- Inspectez l'ensemble des raccordements fluidiques sur la tête de pompe afin d'y détecter d'éventuels signes de fuite. Lorsque la pression du système est redescendue à zéro, resserrez les raccords non étanches.
- Vérifiez s'il y a des fuites dans le système de rinçage des joints. Reportez-vous à la section Test d'étanchéité du système de rinçage des joints () page 282).
- 13. En fonction de la situation :

Situation	Description
Réinstallation de la tête de pompe	Vérifiez s'il y a des fuites dans la pompe. Exécutez le script : E04 – Test de fuite détaillé (pompe) .
Installation d'une nouvelle tête de pompe (tête de pompe de remplacement)	 Réglez les capteurs de pression. Exécutez le script : C05 – Régler les capteurs de pression du système. Vérifiez s'il y a des fuites dans la pompe. Exécutez le script : E04 – Test de fuite détaillé (pompe).

Consultez également

- Retrait de la tête de pompe (> page 197)
- Purge (vue d'ensemble) (▶ page 65)

7.5.3 Remplacement des pistons

À quel moment exécuter cette opération ?

Piston endommagé

Procédez comme suit

- Retirez le piston. Reportez-vous à la section Retrait des pistons (▶ page 201).
- Installez le piston. Reportez-vous à la section Installation des pistons (> page 202).

7.5.3.1 Retrait des pistons

À quel moment exécuter cette opération ?

- Pour remplacer un piston
- Pour effectuer des travaux de maintenance sur les éléments de la tête de pompe

Préparatifs

Retirez la tête de pompe. Reportez-vous à la section Retrait de la tête de pompe (> page 197).

Procédez comme suit

Avec vos doigts, sortez le piston que vous souhaitez retirer de la tête de pompe. Si vous retirez le piston pendant la maintenance des éléments de la tête de pompe, attrapez le piston par sa base et évitez de toucher la surface du piston.



Illustration 92: Retrait des pistons

Consultez également

Installation des pistons (▶ page 202)

7.5.3.2 Installation des pistons

À quel moment exécuter cette opération ?

- Pour installer un piston de remplacement après le retrait du piston
- Pendant la maintenance des éléments de la tête de pompe

Pièces et éléments complémentaires nécessaires

- Piston
- Isopropanol

Préparatifs

Lorsque vous réinstallez un piston que vous avez retirez, nettoyez-le. Reportez-vous à la section Nettoyage du piston (> page 203).

Procédez comme suit



Illustration 93: Installation des pistons

N°	Description
1	Cavité du piston
2	Base du piston

- À l'aide d'une pipette, injectez quelques gouttes d'isopropanol dans la cavité de piston dans laquelle vous souhaitez installer le piston. Lorsque vous réinstallez un piston, installez le piston dans la position où il était lors du retrait.
- 2. Attrapez le piston par sa base.
- 3. Insérez le piston aussi loin que possible dans la cavité de piston.
- Installez la tête de pompe. Reportez-vous à la section Installation de la tête de pompe (▶ page 198).

Consultez également

■ Retrait des pistons (▶ page 201)

7.5.4 Nettoyage du piston

À quel moment exécuter cette opération ?

Avant de réinstaller un piston que vous avez retiré

Matériel nécessaire

- Serviette en papier sèche, non pelucheuse
- Isopropanol

Procédez comme suit

- 1. Inspectez le piston pour vérifier qu'il n'est pas endommagé.
- 2. Rincez soigneusement le piston avec de l'isopropanol.
- 3. À l'aide d'une serviette en papier sèche et non pelucheuse, frottez le piston plusieurs fois.
- 4. Maintenez le piston et passez doucement votre ongle sur la surface du piston. Vous ne devez pas trouver de zones rugueuses ou inégales.
- 5. En fonction de la situation :

Situation	Action
Aucun dommage visible Aucune zone rugueuse ou inégale	 Rincez de nouveau soigneusement le piston avec de l'isopropanol. À l'aide d'une serviette en papier sèche et non pelucheuse, frottez le piston plusieurs fois.
	3. Réinstallez le piston.
Le piston est rayé ou strié	Installez un nouveau piston.

Consultez également

Installation des pistons (▶ page 202)

7.5.5 Remplacement de la bague de support

À quel moment exécuter cette opération ?

Dommages sur la bague de support

Procédez comme suit

- 1. Retirez la bague de support. Reportez-vous à la section Retrait de la bague de support (▶ page 204).
- 2. Installez la bague de support. Reportez-vous à la section Installation de la bague de support (▶ page 206).

7.5.5.1 Retrait de la bague de support

À quel moment exécuter cette opération ?

Pour remplacer la bague de support

Outils nécessaires

- Tournevis type Torx T30
- Outil de manipulation des joints



Illustration 94: Outil de manipulation des joints (à gauche : outil fermé ; à droite : outil ouvert)

N°	Description
1	Broche de retrait
2	Broche d'insertion
3	Molette pour l'ouverture et la fermeture de l'outil de manipulation des joints (l'ouverture et la fermeture sont requises seulement en utilisant la broche de retrait)

Préparatifs

- Retirez la tête de pompe. Reportez-vous à la section Retrait de la tête de pompe (▶ page 197).
- Retirez le piston dont vous souhaitez remplacer la baque de support. Reportez-vous à la section Retrait des pistons (▶ page 201).

Éléments touchés sur la tête de pompe



Illustration 95: Bague de support (ici : pour le piston gauche)

N°	Description
1	Tête de pompe
2	Bague de support
3	Joints de la tête de pompe, (pour la plaque de rinçage des joints et ses vis)
4	Plaque de rinçage des joints
5	Vis, plaque de rinçage des joints (4 vis)
6	Piston

Procédez comme suit

1. À l'aide du tournevis, retirez les vis sur la plaque de rinçage des joints.



Illustration 96: Vis sur la plaque de rinçage des joints (ici : plaque gauche)

- 2. Retirez la plaque de rinçage des joints de la tête de pompe.
- 3. Si un ou plusieurs joints se détachent quand vous retirez la plaque de rinçage des joints, gardez-les pour les réinstaller.

4. Retirez la bague de support : Enfoncez de manière légèrement inclinée la broche *d'insertion* de l'outil de manipulation des joints dans la tête de pompe. N'utilisez pas la broche de *retrait*.



Illustration 97: Bague de support (ici : pour le piston gauche)

5. Retirez l'outil. La bague de support est ainsi enlevée en même temps que l'outil.

Consultez également

Installation de la bague de support (▶ page 206)

7.5.5.2 Installation de la bague de support

À quel moment exécuter cette opération ?

Pour installer une bague de support de remplacement après le retrait de la bague de support

Éléments nécessaires

- Bague de support
- Joints de la tête de pompe pour réinstallation, le cas échéant

Outils nécessaires

- Tournevis type Torx T30
- Outil de manipulation des joints



Illustration 98: Outil de manipulation des joints (à gauche : outil fermé ; à droite : outil ouvert)

N°	Description
1	Broche de retrait
2	Broche d'insertion
3	Molette pour l'ouverture et la fermeture de l'outil de manipulation des joints (l'ouverture et la fermeture sont requises seulement en utilisant la broche de retrait)

Éléments touchés sur la tête de pompe



Illustration 99: Bague de support (ici : pour le piston gauche)

N°	Description
1	Tête de pompe
2	Bague de support
3	Joints de la tête de pompe, (pour la plaque de rinçage des joints et ses vis)
4	Plaque de rinçage des joints
5	Vis, plaque de rinçage des joints (4 vis)
6	Piston

Procédez comme suit

1. Glissez la bague de support sur la broche *d'insertion* de l'outil de manipulation des joints.



Illustration 100: Bague de support glissée sur la broche d'insertion

N°	Description
1	Bague de support
2	Broche d'insertion, outil de manipulation des joints

2. À l'aide de la broche d'insertion, enfoncez de l'outil de manipulation des joints dans la tête de pompe puis retirez l'outil. La bague de support reste dans la tête de pompe.



Illustration 101: Bague de support installée (ici : pour le piston gauche)

- Vérifiez que les joints de la tête de pompe sont correctement installés. Si les joints se sont détachés ou s'ils ne sont pas bien installés dans les fentes de réception, réinstallez-les. Reportez-vous à la section Installation des joints de la tête de pompe (▶ page 222).
- 4. Placez la plaque de rinçage des joints sur la tête de pompe et insérez les vis.



Illustration 102: Vis sur la plaque de rinçage des joints (ici : plaque gauche)

- 5. À l'aide du tournevis, serrez les vis en diagonale.
- 6. Serrez une nouvelle fois les vis en diagonale.
- Installez le piston. Reportez-vous à la section Installation des pistons (▶ page 202).
- Installez la tête de pompe. Reportez-vous à la section Installation de la tête de pompe (▶ page 198).

Consultez également

Retrait de la bague de support (> page 204)

7.5.6 Remplacement du joint du piston

À quel moment exécuter cette opération ?

- Joint de piston endommagé, joint qui fuit
- Débits instables ou bruit de ligne de base

Procédez comme suit

- 1. Retirez le joint de piston. Reportez-vous à la section Retrait du joint du piston (▶ page 209).
- 2. Installez le joint du piston. Reportez-vous à la section Installation du joint de piston (▶ page 212).

7.5.6.1 Retrait du joint du piston

À quel moment exécuter cette opération ?

Pour remplacer le joint du piston

Outils nécessaires

- Tournevis type Torx T30
- Outil de manipulation des joints



Illustration 103: Outil de manipulation des joints (à gauche : outil fermé ; à droite : outil ouvert)

N°	Description
1	Broche de retrait
2	Broche d'insertion
3	Molette pour l'ouverture et la fermeture de l'outil de manipulation des joints (l'ouverture et la fermeture sont requises seulement en utilisant la broche de retrait)

Préparatifs

- Retirez la tête de pompe. Reportez-vous à la section Retrait de la tête de pompe (> page 197).
- 2. Retirez le piston dont vous souhaitez remplacer le joint. Reportezvous à la section Retrait des pistons (▶ page 201).

Éléments touchés sur la tête de pompe



Illustration 104: Joint du piston (ici : pour le piston gauche)

N°	Description
1	Tête de pompe
2	Joint de piston
3	Bague de support
4	Joints de la tête de pompe, (pour la plaque de rinçage des joints et ses vis)
5	Plaque de rinçage des joints
6	Vis, plaque de rinçage des joints (4 vis)
7	Piston

Procédez comme suit

1. À l'aide du tournevis, retirez les vis sur la plaque de rinçage des joints.



Illustration 105: Vis sur la plaque de rinçage des joints (ici : plaque gauche)

- 2. Retirez la plaque de rinçage des joints de la tête de pompe.
- 3. Si un ou plusieurs joints se détachent quand vous retirez la plaque de rinçage des joints, gardez-les pour les réinstaller.

4. Retirez la bague de support : Enfoncez de manière légèrement inclinée la broche *d'insertion* de l'outil de manipulation des joints dans la tête de pompe. N'utilisez pas la broche de *retrait*.



Illustration 106: Bague de support (ici : pour le piston gauche)

5. Retirez l'outil. La bague de support est ainsi enlevée en même temps que l'outil.

Gardez la bague de support pour la réinstaller. Elle ne fait pas partie des pièces d'usure et elle n'a pas besoin d'être remplacée quand vous remplacez le joint de piston.

6. Ouvrez l'outil de manipulation des joints et enfoncez la broche de *retrait* aussi loin que possible dans la tête de pompe.



Illustration 107: Joint du piston (ici : pour le piston gauche)

- 7. Fermez l'outil de manipulation des joints et retirez-le de la tête de pompe. Le joint est ainsi enlevé en même temps que l'outil.
- 8. Pour retirer le joint de l'outil de manipulation des joints : ouvrez l'outil et enlevez le joint de l'outil. Une fois le joint retiré, il ne peut plus être utilisé.

Consultez également

Installation du joint de piston (▶ page 212)

7.5.6.2 Installation du joint de piston

À quel moment exécuter cette opération ?

Pour installer un joint de piston de remplacement après le retrait du joint

Éléments nécessaires

- Joint de piston
- Bague de support retirée pendant le retrait du joint de piston La bague de support ne fait pas partie des pièces d'usure et peut donc être réinstallée. Installez une nouvelle bague de support seulement quand elle est endommagée.
- Joints de la tête de pompe pour réinstallation, le cas échéant

Outils et éléments complémentaires nécessaires

- Tournevis type Torx T30
- Seringue remplie d'isopropanol
- Écouvillon de nettoyage
- Serviette en papier non pelucheuse
- Outil de manipulation des joints



Illustration 108: Outil de manipulation des joints (à gauche : outil fermé ; à droite : outil ouvert)

N°	Description
1	Broche de retrait
2	Broche d'insertion
3	Molette pour l'ouverture et la fermeture de l'outil de manipulation des joints (l'ouverture et la fermeture sont requises seulement en utilisant la broche de retrait)

Éléments touchés sur la tête de pompe



Illustration 109: Joint du piston (ici : pour le piston gauche)

N°	Description
1	Tête de pompe
2	Joint de piston
3	Bague de support
4	Joints de la tête de pompe, (pour la plaque de rinçage des joints et ses vis)
5	Plaque de rinçage des joints
6	Vis, plaque de rinçage des joints (4 vis)
7	Piston

Procédez comme suit

- 1. Nettoyez la bague de support à l'intérieur et à l'extérieur, à l'aide d'isopropanol, d'un écouvillon de nettoyage et d'une serviette en papier non pelucheuse.
- 2. À l'aide d'une pipette, déposez quelques gouttes d'isopropanol dans la cavité de la tête de pompe, sur le bord devant recevoir le joint de piston.



Illustration 110: Bord devant recevoir le joint de piston

3. Glissez la bague de support et le joint de piston sur la broche d'insertion de l'outil de manipulation des joints. Le côté ouvert du joint de piston est orienté vers l'extérieur de l'outil.



Illustration 111: Bague de support et joint de piston glissés sur la broche d'insertion

N°	Description
1	Joint de piston (côté ouvert orienté vers l'extérieur de l'outil)
2	Bague de support
3	Broche d'insertion, outil de manipulation des joints

4. À l'aide de la broche, enfoncez de l'outil de manipulation des joints dans la tête de pompe puis retirez l'outil. Le joint de piston et la bague de support restent dans la tête de pompe.



Illustration 112: Bague de support installée, couvrant le joint de piston (ici : pour le piston gauche)

 Vérifiez que les joints de la tête de pompe sont correctement installés. Si les joints se sont détachés ou s'ils ne sont pas bien installés dans les fentes de réception, réinstallez-les. Reportez-vous à la section Installation des joints de la tête de pompe (▶ page 222). 6. Placez la plaque de rinçage des joints sur la tête de pompe et insérez les vis.



Illustration 113: Vis sur la plaque de rinçage des joints (ici : plaque gauche)

- 7. À l'aide du tournevis, serrez les vis en diagonale.
- 8. Serrez une nouvelle fois les vis en diagonale.
- 9. Installez le piston. Reportez-vous à la section Installation des pistons (> page 202).
- 10. Installez la tête de pompe. Reportez-vous à la section Installation de la tête de pompe (> page 198). Cependant, avant d'exécuter le script pour tester la pompe à la recherche de fuites, faites fonctionner la pompe pendant 30 minutes avec une contre-pression de 80 MPa (recommandé) pour permettre aux nouveaux joints de piston de fonctionner.

Consultez également

Retrait du joint du piston (> page 209)

7.5.7 Remplacement du joint de rinçage des joints

À quel moment exécuter cette opération ?

Joint de rinçage des joints endommagé, observation de liquide en dessous de la tête de pompe (joint qui fuit)

Procédez comme suit

- Retirez le joint de rinçage des joints. Reportez-vous à la section Retrait du joint de rinçage des joints (▶ page 216).
- 2. Installez le joint de rinçage des joints. Reportez-vous à la section Installation du joint de rinçage des joints (▶ page 218).

7.5.7.1 Retrait du joint de rinçage des joints

À quel moment exécuter cette opération ?

Pour remplacer le joint de rinçage des joints

Outils nécessaires

Outil de manipulation des joints



Illustration 114: Outil de manipulation des joints (à gauche : outil fermé ; à droite : outil ouvert)

N°	Description
1	Broche de retrait
2	Broche d'insertion
3	Molette pour l'ouverture et la fermeture de l'outil de manipulation des joints (l'ouverture et la fermeture sont requises seulement en utilisant la broche de retrait)

Préparatifs

- Retirez la tête de pompe. Reportez-vous à la section Retrait de la tête de pompe (> page 197).
- Retirez le piston dont vous souhaitez remplacer le joint de rinçage des joints. Reportez-vous à la section Retrait des pistons (> page 201).
Éléments touchés sur la tête de pompe



Illustration 115: Joint de rinçage des joints (ici : pour le piston gauche)

N°	Description
1	Tête de pompe
2	Plaque de rinçage des joints
3	Joint de rinçage des joints
4	Piston

Procédez comme suit

1. Ouvrez l'outil de manipulation des joints et enfoncez la broche de *retrait* aussi loin que possible dans la plaque de rinçage des joints.



Illustration 116: Joint de rinçage des joints (ici : pour le piston gauche)

- Fermez l'outil de manipulation des joints et retirez-le de la plaque de rinçage des joints. Le joint est ainsi enlevé en même temps que l'outil.
- 3. Pour retirer le joint de l'outil de manipulation des joints : ouvrez l'outil et enlevez le joint de l'outil. Une fois le joint retiré, il ne peut plus être utilisé.

Consultez également

Installation du joint de rinçage des joints () page 218)

7.5.7.2 Installation du joint de rinçage des joints

À quel moment exécuter cette opération ?

Pour installer un joint de rinçage des joints de remplacement après le retrait du joint

Pièces et outils nécessaires

- Joint de rinçage des joints
- Outil de manipulation des joints



Illustration 117: Outil de manipulation des joints (à gauche : outil fermé ; à droite : outil ouvert)

N°	Description
1	Broche de retrait
2	Broche d'insertion
3	Molette pour l'ouverture et la fermeture de l'outil de manipulation des joints (l'ouverture et la fermeture sont requises seulement en utilisant la broche de retrait)

Éléments touchés sur la tête de pompe



Illustration 118: Joint de rinçage des joints (ici : pour le piston gauche)

N°	Description
1	Tête de pompe
2	Plaque de rinçage des joints
3	Joint de rinçage des joints
4	Piston

1. Glissez le joint de rinçage des joints sur la broche d'insertion de l'outil de manipulation des joints. Notez bien l'orientation du joint.



Illustration 119: Joint de rinçage des joints glissant sur la broche d'insertion

N°	Description
1	Joint de rinçage des joints (la flèche indique le côté qui doit être tourné vers l'outil)
2	Broche d'insertion, outil de manipulation des joints

2. À l'aide de la broche d'insertion, enfoncez de l'outil de manipulation des joints dans la plaque de rinçage des joints puis retirez l'outil. Le joint de rinçage des joints reste dans la plaque de rinçage des joints.



Illustration 120: Joint de rinçage des joints installé (ici : pour le piston gauche)

- Installez le piston. Reportez-vous à la section Installation des pistons (> page 202).
- Installez la tête de pompe. Reportez-vous à la section Installation de la tête de pompe (▶ page 198).

Consultez également

Retrait du joint de rinçage des joints (> page 216)

7.5.8 Remplacement des joints de la tête de pompe

À quel moment exécuter cette opération ?

- Détérioration des joints de tête de pompe
- Oservation de liquide en dessous de la tête de pompe (joint qui fuit)

Procédez comme suit

- Retirez les joints de tête de pompe. Reportez-vous à la section Retrait des joints de la tête de pompe (▶ page 220).
- Installez les joints de tête de pompe. Reportez-vous à la section Installation des joints de la tête de pompe (▶ page 222).

7.5.8.1 Retrait des joints de la tête de pompe

À quel moment exécuter cette opération ?

Pour remplacer les joints de tête de pompe

Outils nécessaires

- Tournevis type Torx T30
- Brucelles

Préparatifs

- Retirez la tête de pompe. Reportez-vous à la section Retrait de la tête de pompe (▶ page 197).
- Retirez les pistons. Reportez-vous à la section Retrait des pistons (▶ page 201).

Éléments touchés sur la tête de pompe



Illustration 121: Joints de la tête de pompe (ici : côté gauche de la tête de pompe)

N°	Description
1	Tête de pompe
2	Joints de tête de pompe, pour les vis de la plaque de rinçage des joints (joints toriques 6,5x1)
3	Joint de la tête de pompe, pour la plaque de rinçage des joints (joints toriques 11x1,5)
4	Plaque de rinçage des joints
5	Vis de la plaque de rinçage des joints
6	Piston

Procédez comme suit

Recommandation : Remplacez tous les joints de tête de pompe en même temps (côté gauche et droit de la tête de pompe).

1. À l'aide du tournevis, retirez les vis sur la plaque de rinçage des joints.



Illustration 122: Vis sur la plaque de rinçage des joints (ici : plaque gauche)

2. Retirez la plaque de rinçage des joints de la tête de pompe.

3. Retirez tous les joints de tête de pompe. L'utilisation de brucelles facilite la procédure.



Illustration 123: Joints de la tête de pompe (ici : côté gauche de la tête de pompe)

N°	Description
1	Joints de la tête de pompe, pour les vis de la plaque de rinçage des joints (joints toriques 6,5x1), restant dans la tête de pompe
2	Joint de la tête de pompe, pour la plaque de rinçage des joints (joints toriques 11x1,5), restant dans la plaque de rinçage des joints, le côté vers la tête de pompe

Consultez également

Installation des joints de la tête de pompe (> page 222)

7.5.8.2 Installation des joints de la tête de pompe

À quel moment exécuter cette opération ?

- Pour installer des joints de tête de pompe de remplacement après le retrait du joint
- Pour réinstaller un joint de la tête de pompe qui s'est détaché pendant la maintenance des éléments de la tête de pompe

Pièces et outils nécessaires

- Joints de tête de pompe
- Tournevis type Torx T30
- Seringue remplie d'isopropanol
- Serviette en papier non pelucheuse

Éléments touchés sur la tête de pompe



Illustration 124: Joints de la tête de pompe (ici : côté gauche de la tête de pompe)

N°	Description
1	Tête de pompe
2	Joints de tête de pompe, pour les vis de la plaque de rinçage des joints (joints toriques 6,5x1)
3	Joint de la tête de pompe, pour la plaque de rinçage des joints (joints toriques 11x1,5)
4	Plaque de rinçage des joints
5	Vis de la plaque de rinçage des joints
6	Piston

Procédez comme suit

- 1. Seulement lors de la réinstallation d'un joint de la tête de pompe : nettoyez le joint de la tête de pompe à l'aide d'isopropanol et d'une serviette en papier non pelucheuse.
- 2. Place le joint de la tête de pompe dans la rainure réceptrice.



Illustration 125: Rainures réceptrices des joints de la tête de pompe (ici : côté gauche de la tête de pompe), dans la tête de pompe et la plaque de rinçage des joints

N°	Description
1	Rainures réceptrices des joints de la tête de pompe dans la tête de pompe ; joints pour les vis de la plaque de rinçage des joints (joint torique 6,5x1)
2	Rainure réceptrice du joint de la tête de pompe dans la plaque de rinçage des joints (côté vers la tête de pompe) ; joint pour la plaque de rinçage des joints (joint torique 11x1,5)

3. Avec vos doigts, enfoncez le joint de la tête de pompe aussi loin que possible dans la rainure.



Illustration 126: Joints de la tête de pompe installés, dans la tête de pompe et la plaque de rinçage des joints (ici : côté gauche de la tête de pompe)

4. Placez la plaque de rinçage des joints sur la tête de pompe et insérez les vis.



Illustration 127: Vis sur la plaque de rinçage des joints (ici : plaque gauche)

- 5. À l'aide du tournevis, serrez les vis en diagonale.
- 6. Serrez une nouvelle fois les vis en diagonale.
- 7. Installez le piston. Reportez-vous à la section Installation des pistons (> page 202).
- 8. Installez la tête de pompe. Reportez-vous à la section Installation de la tête de pompe (▶ page 198).

Consultez également

Retrait des joints de la tête de pompe () page 220)

7.5.9 Entretien des clapets de non-retour de la tête de pompe

La tête de pompe est munie de deux clapets de non-retour : un clapet de non-retour d'aspiration et un clapet de non-retour de refoulement.



Illustration 128: Clapets de non-retour de la tête de pompe

N°	Description
1	Clapet de non-retour d'aspiration
2	Clapet de non-retour de refoulement

Suivez les instructions de la procédure de maintenance que vous souhaitez effectuer.

- Pour remplacer le clapet de non-retour d'aspiration, reportez-vous à la section Remplacement du clapet de non-retour d'aspiration (tête de pompe) (> page 226).
- Pour remplacer le clapet de non-retour de refoulement, reportezvous à la section Remplacement du clapet de non-retour de refoulement (tête de pompe) (▶ page 228).
- Pour nettoyer les clapets de non-retour, reportez-vous à la section Nettoyage des clapets de non-retour (tête de pompe) (> page 229).

NOTE Un script qui vous guide pour la maintenance des clapets de nonretour est également disponible. Exécutez le script : **C10 – Nettoyer ou remplacer les clapets de non-retour de la tête de pompe**.

7.5.9.1 Remplacement du clapet de non-retour d'aspiration (tête de pompe)

À quel moment exécuter cette opération ?

- Problème avec le clapet de non-retour ne pouvant pas être résolu par un nettoyage (par exemple, une fuite persistante)
- Clapet de non-retour d'aspiration endommagé

Éléments nécessaires

Clapet de non-retour d'aspiration

Outils et éléments complémentaires nécessaires

- Solvant convenant à la purge, afin d'évacuer toute substance nocive
- Solvant convenant à votre application

Éléments touchés sur la tête de pompe



Illustration 129: Clapet de non-retour d'aspiration sur la tête de pompe

N°	Description
1	Conduite de solvant
2	Clapet de non-retour d'aspiration
3	Tête de pompe

Préparatifs

1. Afin d'évacuer toutes les substances nocives, purgez la pompe à l'aide d'un solvant approprié.

Procédez comme suit

- 1. Sur le clapet de non-retour d'aspiration, débranchez la conduite de solvant.
- 2. Desserrez et retirez le clapet de non-retour d'aspiration avec vos doigts.
- 3. Resserrez le nouveau clapet de non-retour d'aspiration avec vos doigts.

- 4. Raccordez la conduite de solvant sur le clapet de non-retour d'aspiration.
- 5. Purgez la pompe à l'aide du solvant de votre application.
- Inspectez le clapet de non-retour et les raccordements afin d'y détecter une éventuelle fuite. Resserrez les raccords non étanches seulement quand il n'y a plus de pression dans le système.
- Vérifiez s'il y a des fuites dans la pompe. Exécutez le script : E04 Test de fuite détaillé (pompe).

Consultez également

Purge (vue d'ensemble) (▶ page 65)

7.5.9.2 Remplacement du clapet de non-retour de refoulement (tête de pompe)

À quel moment exécuter cette opération ?

- Problème avec le clapet de non-retour ne pouvant pas être résolu par un nettoyage (par exemple, une fuite persistante)
- Clapet de non-retour de refoulement endommagé

Éléments nécessaires

Clapet de non-retour de refoulement

Outils et éléments complémentaires nécessaires

- Solvant convenant à la purge, afin d'évacuer toute substance nocive
- Solvant convenant à votre application

Éléments touchés sur la tête de pompe



Illustration 130: Clapet de non-retour de refoulement sur la tête de pompe

N°	Description
1	Tête de pompe

N°	Description
2	Clapet de non-retour de refoulement
3	Tube en U

Préparatifs

1. Afin d'évacuer toutes les substances nocives, purgez la pompe à l'aide d'un solvant approprié.

Procédez comme suit

- 1. Sur la tête de pompe, retirez le tube en U.
- 2. Desserrez et retirez le clapet de non-retour de refoulement avec vos doigts.
- 3. Insérez et resserrez le nouveau clapet de non-retour de refoulement avec vos doigts.
- 4. Raccordez le tube en U sur la tête de pompe.
- 5. Purgez la pompe à l'aide du solvant convenant à votre application.
- Inspectez le clapet de non-retour et les raccordements afin d'y détecter une éventuelle fuite. Resserrez les raccords non étanches seulement quand il n'y a plus de pression dans le système.
- Vérifiez s'il y a des fuites dans la pompe. Exécutez le script : E04 Test de fuite détaillé (pompe).

Consultez également

Purge (vue d'ensemble) (▶ page 65)

7.5.9.3 Nettoyage des clapets de non-retour (tête de pompe)

La procédure de nettoyage comporte les étapes suivantes :

- Nettoyage à ultrasons avec de l'isopropanol pur
- Rodage des clapets de non-retour de la pompe à une contrepression élevée

À quel moment exécuter cette opération ?

- Mensuellement
- Clapets de non-retour sales ou ne fonctionnant pas correctement

Matériel nécessaire

- Seringue (12 mL), remplie d'isopropanol
- Petit bécher rempli d'isopropanol

- Bain à ultrasons
- Solvants convenant à votre application (pour la purge)
- Solvants convenant à la procédure de rodage
 Dans la plupart des cas, vous pouvez utiliser les solvants de votre
 application. Pour éviter les décharges électrostatiques dans le
 capillaire de contre-pression, utilisez au moins 20 % d'eau dans le
 mélange des solvants A et B à 50/50. Si vous utilisez des solvants
 sans eau comme des solvants de chromatographie en phase
 normale, utilisez 80 % d'isopropanol dans de l'eau dans la voie A et
 dans la voie B.
- Capillaire de contre-pression approprié, générant environ 75 MPa à un débit de 50 μL/min dans l'eau (par exemple, capillaire nanoViper, silice fondue, diamètre intérieur : 20 μm, longueur : 350 mm)

Préparatifs

Retirez les clapets de non-retour que vous souhaitez nettoyer.

Recommandation : Nettoyez les clapets de non-retour d'aspiration et les clapets de non-retour de refoulement de deux têtes de pompe en même temps.

Procédez comme suit

- 1. Dans le sens de l'écoulement, remplissez les clapets de non-retour avec de l'isopropanol à l'aide de la seringue.
- 2. Placez les clapets de non-retour dans le bécher. Nettoyez-les dans le bain à ultrasons pendant 10 minutes.
- 3. Dans la direction du flux, rincez les clapets de non-retour avec de l'isopropanol à l'aide de la seringue.
- 4. Insérez et resserrez les clapets de non-retour. Rebranchez le tube en U ou la conduite de solvant.
- Installez le capillaire de contre-pression sur l'orifice de refoulement de la pompe. Dirigez l'extrémité libre du capillaire vers le bac d'évacuation.
- Laissez la pompe livrer le solvant adapté à la procédure de rodage : Commencez avec 50 % de solvant B et un débit de 10 μL/min. Augmentez le débit jusqu'à ce que la pression du système soit environ 90 MPa et faites fonctionner la pompe pendant 20 minutes.
- Inspectez la tête de pompe et vérifiez si les clapets de non-retour présentent des signes de fuite. Resserrez les raccords non étanches seulement quand il n'y a plus de pression dans le système.

- 8. Désinstallez le capillaire de contre-pression et raccordez le système selon les besoins de votre application.
- 9. Purgez la pompe à l'aide du solvant convenant à votre application.
- 10. Vérifiez s'il y a des fuites dans la pompe. Exécutez le script : **E04 Test de fuite détaillé (pompe)**.

Consultez également

- Remplacement du clapet de non-retour d'aspiration (tête de pompe)
 () page 226)
- Remplacement du clapet de non-retour de refoulement (tête de pompe) (> page 228)

7.6 Maintenance du débitmètre

La présente section décrit comment entretenir ou remplacer les éléments de la tête de pompe. Suivez les instructions de la procédure de maintenance que vous souhaitez effectuer.

7.6.1 Remplacement du filtre du débitmètre

À quel moment exécuter cette opération ?

Filtre obstrué

Le script de diagnostics suggère de remplacer le filtre

Éléments nécessaires

Kit du filtre du débitmètre

Outils et éléments complémentaires nécessaires

- Clé ouverte, taille : 13 mm
- Solvant convenant à l'élimination de substances nocives
- Solvant convenant à votre application

Éléments touchés



Illustration 131: Filtre du débitmètre (à gauche : éléments installés ; à droite : filtre retiré)

N°	Description
1	Tête de pompe
2	Capillaire raccordant la tête de pompe et le débitmètre
3	Filtre du débitmètre

Préparatifs

- 1. Afin d'évacuer toutes les substances nocives, purgez la pompe à l'aide d'un solvant approprié.
- 2. Préparez le filtre de remplacement.

Préparation du filtre de remplacement



Illustration 132: Assemblage du filtre du débitmètre

N°	Description
1	Vis du filtre
2	Joint de la vis du filtre
3	Fritté de filtre

- 1. Avec vos doigts, enfoncez le joint sur la vis du filtre jusqu'à ce qu'il soit placé sur le bord récepteur.
- 2. À l'aide d'une pipette, déposez une goutte de solvant sur la vis du filtre et insérez le fritté du filtre. Le solvant permet de garder le fritté dans la vis.

Procédez comme suit

- 1. Retirez le capillaire qui raccorde la tête de pompe et le débitmètre.
- 2. Avec la clé, desserrez la vis du filtre et retirez le filtre de la pompe.
- 3. Insérez le nouveau filtre et resserrez la vis du filtre avec la clé.
- 4. Branchez le capillaire qui raccorde la tête de pompe et le débitmètre.
- 5. Purgez la pompe et le débitmètre à l'aide d'un solvant convenant à l'application.

- 6. Inspectez les raccordements pour voir s'ils présentent des signes de fuite. Resserrez les raccords non étanches quand il n'y a plus de pression dans le système.
- Vérifiez si le nouveau filtre fonctionne correctement. Exécutez le script : D01 – Tester la contre-pression du système.
- Vérifiez s'il y a des fuites dans la pompe. Exécutez le script : E04 Test de fuite détaillé (pompe).

Consultez également

Purge (vue d'ensemble) (▶ page 65)

7.6.2 Remplacement de la conduite d'évacuation du débitmètre

À quel moment exécuter cette opération ?

Conduite d'évacuation endommagée ou obstruée

Pièces et éléments complémentaires nécessaires

- Conduite d'évacuation du débitmètre
- Solvant convenant à l'élimination de substances nocives

Éléments touchés





N°	Description
1	Conduite d'évacuation du débitmètre
2	Bloc du détecteur de rinçage des joints
3	Orifice d'évacuation
4	Position d'installation de l'assemblage du détecteur de rinçage des joints
5	Orifice de refoulement de l'évacuation (vers le dispositif d'évacuation)

N°	Description
6	Rainure dans la paroi de l'orifice d'évacuation
7	Capteur de fuites

Préparatifs

- 1. Afin d'évacuer toutes les substances nocives, purgez la pompe à l'aide d'un solvant approprié.
- Pour accéder facilement à la paroi de l'orifice d'évacuation et l'orifice de refoulement de l'évacuation, retirez le bloc de détection de rinçage des joints (pied et détecteur).

Retrait de l'assemblage du détecteur de rinçage des joints

Saisissez et tirez le bloc du détecteur de rinçage des joints (pied et détecteur) vers le haut par le pied. Le détecteur incluant le tube et le câble peut rester raccordé au pied.



Illustration 134: Retrait du bloc de détecteur de rinçage des joints

N°	Description
1	Pied, détecteur de rinçage des joints
2	Détecteur de rinçage des joints
3	Câble du détecteur
4	Conduite de rinçage des joints

Procédez comme suit

- 1. Retirez la conduite d'évacuation du port Purge Out du débitmètre et de l'orifice de refoulement de l'évacuation.
- 2. Branchez la nouvelle conduite d'évacuation sur le port Purge Out.

- Dirigez la conduite à travers la rainure dans la paroi de l'orifice d'évacuation dans l'orifice de refoulement de l'évacuation.
 Pour éviter toute fausse alarme de fuite, la conduite d'évacuation doit être bien positionnée dans l'orifice de refoulement de l'évacuation. Sinon, le liquide peut s'accumuler dans l'orifice d'évacuation et activer le capteur de fuite.
- 4. Réinsérez le bloc du détecteur de rinçage des joints.

Consultez également

Purge (vue d'ensemble) (▶ page 65)

7.7 Entretien de la conduite de solvant

La présente section décrit comment entretenir ou remplacer les conduites de solvants ou le filtre de conduite de solvant. Suivez les instructions de la procédure de maintenance que vous souhaitez effectuer.

7.7.1 Vidange des conduites de solvant

À quel moment exécuter cette opération ?

Pour vider les conduites de solvant, par exemple pour des procédures de maintenance ou de service

Matériel nécessaire

Solvant convenant à l'élimination de substances nocives, le cas échéant

Procédez comme suit

- 1. Afin d'évacuer toutes les substances nocives, le cas échéant, purgez la pompe à l'aide d'un solvant approprié.
- 2. Dévissez le bouchon du réservoir de solvant et retirez la conduite de solvant en même temps que le bouchon du réservoir.
- 3. Lancez un cycle de purge, si besoin.
- 4. Arrêtez la purge dès que la conduite de solvant est vide.

Consultez également

Purge (vue d'ensemble) (▶ page 65)

7.7.2 Remplacement des conduites de solvant entre le réservoir et la pompe

À quel moment exécuter cette opération ?

Détérioration ou obstruction de la conduite de solvant

Pièces et éléments complémentaires nécessaires

- Conduites de solvant entre le réservoir de solvant et la pompe
- Solvant convenant à votre application

Éléments touchés



Illustration 135: Conduites de solvant entre le réservoir et la pompe

N°	Description
1	Conduites de solvant des réservoirs, dirigées à la pompe à travers tous les modules situés au-dessus de la pompe dans le système empilé avec
1a	Conduite de solvant, voie A
1b	Conduite de solvant, voie B
2	Vannes d'arrêt, A et B
3	Bloc du détecteur de rinçage des joints

Préparatifs

- 1. Videz les conduites de solvant. Reportez-vous à la section Vidange des conduites de solvant (▶ page 237).
- Si vous voulez remplacer le filtre de conduite de solvant en même temps que la conduite de solvant, préparez le filtre. Reportez-vous à la section Remplacement de filtre de conduite de solvant (> page 243).
- 3. Pour y accéder facilement, retirez le bloc de détection de rinçage des joints (pied et détecteur).

Retrait de l'assemblage du détecteur de rinçage des joints

Saisissez et tirez le bloc du détecteur de rinçage des joints (pied et détecteur) vers le haut par le pied. Le détecteur incluant le tube et le câble peut rester raccordé au pied.





N°	Description
1	Pied, détecteur de rinçage des joints
2	Détecteur de rinçage des joints
3	Câble du détecteur
4	Conduite de rinçage des joints

Retrait de la conduite de solvant

- 1. Retirez la conduite de solvant du bouchon de réservoir :
 - a) Retirez le filtre de la conduite de solvant.
 - b) Retirez le guide de maintien.
 - c) Sortez la conduite de solvant du bouchon de réservoir.
- 2. Débranchez la conduite de solvant dans la pompe.
- Retirez la conduite de solvant de la pompe et tous les modules situés au-dessus de la pompe dans le système empilé. Lorsque vous retirez la conduite de solvant des guides de tubulure, ne tirez pas sur d'autres tubulures présentes dans les guides.

Acheminement de la nouvelle conduite de solvant jusqu'au bac à solvant

1. Branchez chaque conduite de solvant à l'orifice d'admission dans la pompe pour lequel elle est prévue (reportez-vous à l'étiquetage présent sur les conduites de solvant et dans la pompe).

- Acheminez les conduites de solvant de la pompe au bac à solvants, en passant par les guides de tubulure dans la pompe et tous les modules situés au-dessus de la pompe dans le système empilé. Respectez les consignes d'acheminement indiquées à la section Passage des capillaires et des tubes dans le système (▶ page 120).
- Acheminez les conduites de solvant à travers le trou de guidage situé à l'intérieur du bac à solvants et poussez-les dans un ou plusieurs guides de tubulure.





N°	Description
А	Trou de guidage (bac à solvants)
1-4	Guides de tubulure

Raccordement de la conduite de solvant au réservoir de solvant



Illustration 138: Préparation de la conduite de solvant et du bouchon de réservoir

N°	Description
1	Conduite de solvant
2	Guide de maintien
3	Bouchon de réservoir

N°	Description
4	Capuchons protecteurs
5	Filtre de conduite de solvant avec adaptateur de conduite de solvant

- 1. Faites passer chaque conduite de solvant par un guide de maintien.
- 2. Faites passer la conduite de solvant par un trou dans un bouchon pour réservoir de solvant. Le guide de maintien permet d'immobiliser la conduite de solvant dans le réservoir.
- 3. Obturez les trous restants des bouchons de réservoir au moyen de capuchons protecteurs.
- 4. Glissez un filtre de conduite de solvant sur chaque conduite de solvant.
- 5. Serrez le bouchon du réservoir manuellement. Enfoncez le guide de maintien dans le trou du bouchon du réservoir afin d'immobiliser la tubulure dans le bouchon.
- 6. Positionnez les conduites de solvant directement dans les guides de tubulure.
- Contrôlez les conduites de solvant sur l'ensemble du trajet d'écoulement. Assurez-vous que les conduites ne sont, à aucun endroit du trajet d'écoulement, pliées, pincées ou serrées.
- 8. Si applicable, réinsérez le bloc du détecteur de rinçage des joints.
- 9. Purgez la pompe à l'aide du solvant convenant à votre application.

Consultez également

Purge (vue d'ensemble) (▶ page 65)

7.7.3 Remplacement des conduites de solvant entre la vanne d'arrêt et la tête de pompe

À quel moment exécuter cette opération ?

Détérioration ou obstruction de la conduite de solvant

Pièces et éléments complémentaires nécessaires

- Conduites de solvant entre la vanne d'arrêt et la tête de pompe
- Solvant convenant à l'élimination de substances nocives
- Solvant convenant pour votre application

Préparatifs

• Afin d'évacuer toutes les substances nocives, purgez la pompe à l'aide d'un solvant approprié.

Éléments touchés



Illustration 139: Conduites de solvant entre les vannes d'arrêt et les têtes de pompe

N°	Description
1	Vannes d'arrêt, A et B
2	Têtes de pompe, A et B
3	Conduite de solvant entre la vanne d'arrêt B et la tête de pompe B
4	Conduite de solvant entre la vanne d'arrêt A et la tête de pompe A

Procédez comme suit

- 1. Débranchez la conduite de solvant sur la vanne d'arrêt et la tête de pompe.
- 2. Installez la nouvelle conduite de solvant.
- 3. Purgez la pompe à l'aide d'un solvant convenant pour l'application.

Consultez également

Purge (vue d'ensemble) (▶ page 65)

7.7.4 Remplacement de filtre de conduite de solvant

À quel moment exécuter cette opération ?

- Pendant une maintenance régulière
- Perméabilité incomplète du fritté de la conduire de solvant
- Filtre de conduite de solvant endommagé

Pièces et éléments complémentaires nécessaires

- Selon les besoins
 - Fritté de filtre
 - Filtre de conduite de solvant (avec porte-filtres, fritté de filtre et adaptateur de conduite de solvant)
- Solvant convenant à votre application

Éléments touchés



Illustration 140: Filtre de conduite de solvant (à gauche : en remplaçant seulement le fritté de filtre ; à droite : en remplaçant le filtre de conduite de solvant)

N°	Description
1	Adaptateur de conduite de solvant
2	Porte-filtre, partie supérieure
3	Fritté de filtre
4	Porte-filtre, partie inférieure

Préparatifs

- 1. Dévissez le bouchon du réservoir de solvant et retirez la conduite de solvant en même temps que le bouchon du réservoir.
- 2. Retirez le porte-filtre (avec adaptateur) de la conduite de solvant.

Procédez comme suit

Comme requis, suivez les étapes de remplacement du fritté de filtre ou de remplacement du filtre.

Étapes de remplacement du fritté dans le filtre de conduite de solvant

- 1. Ouvrez le porte-filtre et retirez le fritté.
- 2. Montez le filtre de conduite de solvant en revêtant des gants propres et appropriés :
 - a) Placez le fritté dans une position de niveau dans le porte-filtre (partie inférieure).
 - b) Vissez le haut du filtre fermement sur le bas du filtre.
- 3. Glissez le porte-filtre avec adaptateur sur la conduite de solvant.
- 4. Resserrez le bouchon du réservoir manuellement. Le guide de maintien doit rester dans le trou du bouchon de réservoir. Si ce n'est pas le cas, enfoncez le guide de maintien dans le trou afin d'immobiliser la tubulure dans le bouchon.
- 5. Pour éliminer tout air des conduites de solvant, purgez la pompe à l'aide du solvant de l'application.

Étapes de remplacement du filtre de conduite de solvant

- 1. Montez le filtre de conduite de solvant en revêtant des gants propres et appropriés :
 - a) Insérez l'adaptateur de conduite de solvant dans le porte-filtre (partie supérieure).
 - b) Tournez le porte-filtre (partie supérieure) avec adaptateur tête en bas et pressez-le sur une surface appropriée pour que l'adaptateur soit fermement attaché sur le porte-filtre.
 - c) Placez le fritté dans une position de niveau dans le porte-filtre (partie inférieure).
 - d) Vissez le haut du filtre fermement sur le bas du filtre.
- 2. Glissez le porte-filtre avec adaptateur sur la conduite de solvant.
- Resserrez le bouchon du réservoir manuellement. Le guide de maintien doit rester dans le trou du bouchon de réservoir. Si ce n'est pas le cas, enfoncez le guide de maintien dans le trou afin d'immobiliser la tubulure dans le bouchon.
- 4. Purgez la pompe à l'aide du solvant de votre application.

Consultez également

Purge (vue d'ensemble) (▶ page 65)

7.8 Entretien de la boucle d'échantillonnage

À quel moment exécuter cette opération ?

- La boucle d'échantillonnage est obstruée ou endommagée
- Différent volume de boucle d'échantillonnage requis

Éléments nécessaires

Boucle d'échantillonnage de remplacement ou boucle d'échantillonnage d'une autre taille

Procédez comme suit

- 1. Retirez la boucle d'échantillonnage.
- 2. Installez la boucle d'échantillonnage de remplacement.
- Finalisez la procédure de remplacement et sélectionnez la boucle d'échantillonnage correspondante dans le système de gestion de données chromatographiques.

NOTE Un script qui vous donne des instructions pour le remplacement de la boucle d'échantillonnage est également disponible. Si vous avez besoin d'être guidé, exécutez le script : **A07 – Remplacer la boucle d'échantillonnage**.

7.8.1 Composants de la boucle d'échantillonnage

Illustration 141: Unité de boucle d'échantillonnage installée sur l'unité d'aiguille

N°	Description
1	Raccord de boucle d'échantillonnage sur l'unité d'aiguille
2	Boucle d'échantillonnage
3	Plaque de la boucle d'échantillonnage
4	Vis de fixation pour la plaque de la boucle d'échantillonnage
5	Logement dans la plaque de la plate porte-boucle

7.8.2 Retrait de la boucle d'échantillonnage

Outils nécessaires

Tournevis type Torx T10

Préparatifs

- 1. Déplacez l'unité d'aiguille à la position d'entretien.
- 2. Retirez le couvercle isolant.
- 3. Sécurisez l'unité d'aiguille.
- 4. Mettez le passeur d'échantillon hors tension au moyen de son interrupteur d'alimentation principal.

Procédez comme suit

 Desserrez légèrement la vis de fixation à l'aide du tournevis. La vis peut rester légèrement vissée au support de l'unité de boucle d'échantillonnage.

- Déconnectez le raccord de boucle d'échantillonnage de l'unité d'aiguille. Pour un accès plus facile, utilisez le tournevis dynamométrique avec embout Viper.
- 3. Déconnectez le raccord de boucle d'échantillonnage de la vanne d'injection.
- 4. Poussez l'unité de boucle d'échantillonnage légèrement vers le haut, de telle sorte que la vis de fixation corresponde à la partie large de l'ouverture dans la plaque de la boucle d'échantillonnage. Retirez l'unité de boucle d'échantillonnage de l'unité d'aiguille.



Illustration 142: Retrait de la boucle d'échantillonnage

 Si la boucle d'échantillonnage doit être réinstallée : Protégez les raccords ouverts de la boucle d'échantillonnage avec des capuchons.

7.8.3 Installation de la boucle d'échantillonnage

Pièces et outils nécessaires

- Si la boucle d'échantillonnage doit être remplacée : boucle d'échantillonnage de rechange
- Tournevis type Torx T10
- Tournevis dynamométrique avec adaptateur d'extension et embout Viper

Procédez comme suit

- Retirez le capuchon du raccord inférieur de la boucle d'échantillonnage et branchez la boucle d'échantillonnage sur la vanne d'injection. Pour en savoir plus sur les branchements, reportez-vous à la section Attributions des orifices de la vanne du passeur d'échantillon (▶ page 48).
- Retirez le capuchon du raccord supérieur de la boucle d'échantillonnage et branchez la boucle d'échantillonnage sur l'unité d'aiguille.

- 3. Suspendez la plaque de la boucle d'échantillonnage sur le support de l'unité de boucle d'échantillonnage :
 - a) Positionnez l'ouverture de la plaque porte-boucle sur la vis de fixation. Assurez-vous que la vis est desserrée.
 - b) Alignez le logement allongé dans la plaque avec le support.
- Introduisez le raccord de la boucle d'échantillonnage dans l'unité d'aiguille et vissez-le à l'intérieur. Assurez-vous que le capillaire en provenance du raccord de l'unité d'aiguille sorte du raccord en pointant vers 5 heures.
- 5. Serrez le raccorde de la boucle d'échantillonnage sur l'unité d'aiguille à l'aide du tournevis dynamométrique. Après avoir serré, le capillaire devrait pointer vers 6 heures.
- 6. Serrez la vis de fixation de la plaque porte-boucle à l'aide du tournevis.



Illustration 143: Alignement de l'unité de boucle d'échantillonnage sur la vis

7. Guidez le capillaire à la vanne d'injection dans le support de montage du couvercle isolant.



Illustration 144: Capillaire guidé à travers le support de montage du couvercle isolant

 Finalisez la procédure de remplacement. Reportez-vous à la section Finalisation de la procédure de remplacement de la boucle d'échantillonnage (> page 249).

7.8.4 Finalisation de la procédure de remplacement de la boucle d'échantillonnage

Après avoir installé la boucle d'échantillonnage, procédez comme suit :

- 1. Déverrouillez l'unité d'aiguille.
- Vérifiez que la boucle d'échantillonnage se déplace librement quand vous enfoncez l'unité d'aiguille dans le compartiment à échantillons. Si la boucle d'échantillonnage reste bloquée contre le panneau intérieur, retirez-la et réinstallez-la à nouveau en suivant les étapes de cette section. Assurez-vous que la boucle d'échantillonnage est correctement alignée et connecteur avec points capillaire à 6 heures.
- 3. Montez le couvercle isolant.
- Mettez le passeur d'échantillon sous tension au moyen de son interrupteur d'alimentation principal. Le passeur d'échantillon effectue un test automatique. L'aiguille s'initialise et ajuste sa position.
- 5. Au besoin, réglez le paramètre **Nominal Loop Volume** (Volume nominal de la boucle) manuellement pour la boucle d'échantillonnage installée. Respectez les volumes indiqués sur l'étiquette de la boucle d'échantillonnage.

Description	Volume nominal de la boucle	Volume total de la boucle
Boucle d'échantillonnage, 10 μL	10 µL	23 μL
Boucle d'échantillonnage, 25 μL	25 μL	50 μL
Boucle d'échantillonnage, 100 μL	100 μL	130 µL

6. Préparez le passeur d'échantillon pour l'utilisation. Reportez-vous à la section Préparer le système à l'utilisation (▶ page 161).

7.9 Entretien de l'unité d'aiguille

7.9.1 Composant de l'unité d'aiguille et du siège de l'aiguille



Illustration 145: Composants de l'unité d'aiguille et du siège de l'aiguille

N°	Description
1	Vis de verrouillage d'expédition
2	Verrou du pousseur de flacon
3	Plaque de positionnement sur le pousseur de l'unité d'aiguille
4	Vis de fixation du pousseur de flacon
5	Vis de fixation de l'unité d'aiguille sur le chariot de l'aiguille
6	Aiguille
7	Pousseur de flacon
8	Siège de l'aiguille

7.9.2 Verrouillage de l'unité d'aiguille

À quel moment exécuter cette opération ?

Avant une procédure de maintenance ou si l'appareil doit être déplacé

AVIS

Si l'unité d'aiguille n'est pas verrouillée en place pendant les opérations de maintenance ou le déplacement de l'appareil, cela peut détériorer le passeur d'échantillon. Verrouillez l'unité d'aiguille lorsque vous en recevez l'instruction afin d'éviter toute détérioration du passeur d'échantillon.

Outils nécessaires

Tournevis plat (facultatif)

Préparatifs

Retirez le couvercle isolant. Reportez-vous à la section Retrait du couvercle isolant (▶ page 97).

Mise en position d'entretien de l'unité d'aiguille

Appuyez une fois sur le bouton **SERVICE** du panneau de commande pour déplacer l'unité d'aiguille vers le haut depuis le siège de l'aiguille et la mettre dans la première position d'entretien.

Serrage de la vis de verrouillage d'expédition

Au-dessus de l'unité d'aiguille, serrez manuellement la vis de verrouillage d'expédition jusqu'à ce qu'elle ne puisse plus être tournée davantage. L'utilisation d'un tournevis pour écrous à fente peut faciliter la manœuvre.



Illustration 146: Serrage de la vis de verrouillage d'expédition

7.9.3 Déverrouillage de l'unité d'aiguille

Outils nécessaires

Tournevis pour écrous à fente (facultatif)

Desserrage de la vis de verrouillage d'expédition

AVIS

Si la vis de verrouillage d'expédition n'est pas correctement desserrée, le passeur d'échantillon peut ne pas fonctionner correctement ou être endommagé. Assurez-vous toujours que la vis de verrouillage d'expédition est complètement desserrée avant de mettre en marche le passeur d'échantillon.

Desserrez la vis de verrouillage d'expédition qui se trouve au-dessus de l'unité d'aiguille. Avec vos doigts, tournez la vis imperdable dans le sens inverse des aiguilles d'une montre jusqu'à ce qu'elle reste suspendue librement dans le ressort. L'utilisation d'un tournevis plat peut faciliter la manœuvre.



Illustration 147: Desserrage de la vis de verrouillage d'expédition
Vérification de l'unité d'aiguille

Selon les besoins :

État du passeur d'échantillon	Actions
Le passeur d'échantillon a été arrêté au cours de la procédure de maintenance	Vérifiez que l'unité d'aiguille se déplace librement : avancez-la dans le compartiment à échantillons le long de l'entraînement horizontal de l'aiguille. Reportez-vous à l'illustration ci- dessous. Si elle se déplace sans difficulté, alors la vis de verrouillage d'expédition est suffisamment desserrée.
Le passeur d'échantillon est resté allumé pendant la procédure de maintenance	Appuyez sur le bouton SERVICE pour déplacer l'unité d'aiguille dans le compartiment.



Illustration 148: Déplacement de l'unité d'aiguille (passeur d'échantillon hors tension)

Redémarrage de l'unité d'aiguille après sa mise en position d'entretien

Selon les besoins :

État du passeur d'échantillon	Actions
Le passeur d'échantillon a été arrêté au cours de la procédure de maintenance	L'unité d'aiguille est initialisée automatiquement à la mise sous tension du passeur d'échantillon.
Le passeur d'échantillon est resté allumé pendant la procédure de maintenance	Selon la position d'entretien utilisée précédemment :
	 Appuyez sur le bouton SERVICE du panneau de commande et attendez que l'aiguille atteigne la première position d'entretien (aiguille vers le haut).
	 Appuyez à nouveau sur le bouton SERVICE et attendez que l'aiguille atteigne la seconde position d'entretien (aiguille au-dessus de l'orifice de rinçage).
	3. Appuyez une troisième fois sur le bouton SERVICE pour initialiser l'unité d'aiguille après sa mise en position d'entretien. Lorsque l'aiguille se déplace, vérifiez visuellement que la boucle d'échantillonnage peut bouger librement.

7.9.4 Remplacement du siège de l'aiguille et de l'unité de l'aiguille

À quel moment exécuter cette opération ?

En fonction du type de problème, il peut être nécessaire de remplacer le siège de l'aiguille et l'unité d'aiguille ensemble ou séparément.

À quel moment exécuter cette opération ?	Mesure corrective
Le siège de l'aiguille est obstrué	 Exécutez le script C21 – Nettoyer ou remplacer le siège et l'unité d'aiguille. Si le script ne permet pas d'éliminer le blocage : remplacez le siège d'aiguille (reportez-vous à la section Remplacement du siège de l'aiguille (* page 255)).
	3. Vérifiez si l'unité d'aiguille est endommagée.
Le siège de l'aiguille fuit	 Serrez le siège de l'aiguille avec un tournevis dynamométrique avec embout pour le siège de l'aiguille. Reportez-vous également à la section Inspection des vannes du passeur d'échantillon à la recherche de fuites (> page 301).
	Si le siège de l'aiguille continue de fuir : Remplacez le siège de l'aiguille et l'unité d'aiguille ensemble.
	 Remplacez le siège de l'aiguille, reportez-vous à la section Remplacement du siège de l'aiguille () page 255).
	3. Remplacez l'unité d'aiguille, reportez-vous à la section Remplacement de l'unité d'aiguille (▶ page 258).

À quel moment exécuter cette opération ?	Mesure corrective
L'aiguille est endommagée	Remplacez le siège de l'aiguille et l'unité d'aiguille ensemble : 1. Remplacez le siège de l'aiguille, reportez-vous à la section Remplacement du siège de l'aiguille (▶ page 255).
	 Remplacez l'unité d'aiguille, reportez-vous à la section Remplacement de l'unité d'aiguille (page 258).

Éléments nécessaires

- Unité d'aiguille
- Siège de l'aiguille

Outils nécessaires

- Tournevis type Torx T10
- Tournevis dynamométrique avec embout pour le siège de l'aiguille



Illustration 149: Outil pour resserrer le siège de l'aiguille

N°	Description
1	Embout pour le siège de l'aiguille
2	Base du tournevis dynamométrique

NOTE Vous pouvez exécuter un script afin d'obtenir des instructions étape par étape directement sur l'appareil. Pour plus d'informations, exécutez le script : **C21 – Nettoyer ou remplacer le siège et l'unité d'aiguille.**

7.9.4.1 Remplacement du siège de l'aiguille

Procédez comme suit

1. Appuyez deux fois sur le bouton **SERVICE** pour mettre l'unité d'aiguille dans la seconde position d'entretien au-dessus de l'orifice de rinçage.

- 2. Mettez le passeur d'échantillon hors tension au moyen de son interrupteur d'alimentation principal.
- 3. Retirez le couvercle isolant.
- 4. Retirez le capillaire du support de montage du couvercle isolant.
- 5. Pincez le support de montage du couvercle isolant et poussez-le vers le haut pour le retirer.



Illustration 150: Retrait du support de montage du couvercle isolant

AVIS—Le tournevis dynamométrique avec l'embout pour le siège de l'aiguille peut être endommagé s'il n'est pas utilisé correctement.

- Insérez l'embout pour le siège de l'aiguille complètement dans le tournevis dynamométrique.
- Insérez le tournevis dynamométrique avec l'embout pour le siège de l'aiguille droit dans le siège de l'aiguille.
- Ne poussez et ne tirez pas sur l'outil avec une force inutile.
- 6. Insérez l'embout pour le siège de l'aiguille dans le tournevis dynamométrique.
- 7. Positionnez le tournevis dynamométrique avec l'embout pour le siège de l'aiguille dans le siège de l'aiguille.

8. Tournez l'outil dans le sens contraire des aiguilles d'une montre pour retirer le siège de l'aiguille.



Illustration 151: Retrait du siège de l'aiguille avec le tournevis dynamométrique

- 9. Déballez le siège d'aiguille de remplacement.
- 10. Insérez le siège de l'aiguille et tournez-le manuellement dans le sens des aiguilles d'une montre.



Illustration 152: Insertion du siège de l'aiguille

 Utilisez le tournevis dynamométrique avec l'embout pour le siège de l'aiguille et tournez dans le sens des aiguilles d'une montre pour serrer le siège de l'aiguille jusqu'à la limite de couple, signalée par un clic.

AVIS—Le siège de l'aiguille pourrait être endommagé par l'utilisation d'autres outils

Le tournevis dynamométrique avec l'embout pour le siège de l'aiguille est destiné au serrage du siège de l'aiguille. À l'aide de cet outil, le siège de l'aiguille est correctement serré sans serrage excessif. N'utilisez pas un autre outil.



Illustration 153: Serrage du siège de l'aiguille

 En fonction du problème, suivez les consignes données aux sections Remplacement de l'unité d'aiguille (▶ page 258) ou Finalisation de la procédure de remplacement du siège de l'aiguille et de l'unité d'aiguille (▶ page 262).

7.9.4.2 Remplacement de l'unité d'aiguille

NOTE Toute détérioration de l'unité d'aiguille se répercute sur le siège de l'aiguille et le détériore aussi. Pour cette raison, il est recommandé de remplacer le siège de l'aiguille et l'unité d'aiguille en même temps.

Procédez comme suit

- Retirez la boucle d'échantillonnage. Suivez les étapes de retrait correspondantes à la section Retrait de la boucle d'échantillonnage (> page 246). Le raccord de la vanne d'injection peut rester en place.
- 2. Déplacez manuellement l'unité d'aiguille un peu vers l'avant.

- 3. Sécurisez l'unité d'aiguille.
- 4. À l'aide du tournevis, desserrez les deux vis qui fixent l'unité d'aiguille sur le chariot de l'aiguille.



Illustration 154: Dévissage de l'unité d'aiguille

N°	Description
1	Vis de fixation de l'unité d'aiguille sur le chariot de l'aiguille

5. Tirez l'unité d'aiguille légèrement vers l'avant et retirez-la du chariot de l'aiguille.

AVIS—L'ouverture de la plaque de positionnement est sensible à la poussière et aux débris.

Veillez à éviter toute pénétration de poussière ou de débris à l'intérieur du passeur d'échantillon.



AVERTISSEMENT—Aiguille d'injection à pointe acérée

L'aiguille d'injection est dotée d'une pointe très acérée qui peut causer des blessures cutanées.

Pour éviter les blessures, ne touchez jamais la pointe de l'aiguille.

AVIS—L'aiguille dans l'unité d'aiguille n'est pas fixée.

- Retirez doucement l'unité d'aiguille après avoir dévissé les vis.
- Manipulez l'unité d'aiguille de remplacement avec soin, en particulier après avoir retiré le verrouillage d'expédition.
- 6. Déballez l'unité d'aiguille de remplacement.
- Utilisez l'unité d'aiguille avec le verrouillage d'expédition monté. Le verrouillage d'expédition se débloquera automatiquement pendant le processus de montage de l'unité d'aiguille et il peut être retiré par la suite.
- 8. Positionnez l'unité d'aiguille de remplacement sur le chariot de l'aiguille. Respectez les consignes suivantes :
 - Le rebord qui dépasse à l'arrière de l'unité d'aiguille doit s'insérer correctement dans l'ouverture du chariot de l'aiguille (reportez-vous à l'illustration).
 - La partie en saillie sur la face arrière de l'unité d'aiguille doit s'ajuster dans le renfoncement du chariot de l'aiguille.



Illustration 155: Installation de l'unité d'aiguille sur le chariot de l'aiguille

N°	Description
1	Ouverture du chariot de l'aiguille
2	Rebord en saillie sur l'unité d'aiguille

9. Placez et serrez les deux vis situées sur l'unité d'aiguille.

- 10. Vérifiez si le verrou et le pousseur de flacon se déplacent facilement vers le haut.



N°	Description
1	Verrou du pousseur de flacon
2	Pousseur de flacon

- 11. Déverrouillez l'unité d'aiguille.
- 12. Déplacez l'unité d'aiguille un peu vers l'arrière au-dessus de l'orifice de rinçage.
- Suivez les consignes indiquées à la section Finalisation de la procédure de remplacement du siège de l'aiguille et de l'unité d'aiguille (▶ page 262).

7.9.4.3 Finalisation de la procédure de remplacement du siège de l'aiguille et de l'unité d'aiguille

1. Réinstallez le support de montage du couvercle isolant en le poussant dans les rails de guidage jusqu'à ce qu'il s'enclenche en position.



Illustration 157: Installation du support de montage du couvercle isolant

- 2. Déplacez manuellement l'unité d'aiguille un peu vers l'avant.
- 3. Sécurisez l'unité d'aiguille.
- Si la boucle d'échantillonnage a été retirée, réinstallez-la. Suivez les étapes d'installation correspondantes figurant à la section Installation de la boucle d'échantillonnage (▶ page 247).
- 5. Guidez le capillaire depuis la boucle d'échantillonnage vers la vanne d'injection à travers le support de montage du couvercle isolant.
- 6. Déverrouillez l'unité d'aiguille.
- 7. Remontez le couvercle isolant.
- 8. Remettez le passeur d'échantillon sous tension au moyen de son interrupteur d'alimentation principal. Le passeur d'échantillon effectue un test automatique. L'aiguille s'initialise et ajuste sa position.
- En fonction du remplacement effectué : Réglez les paramètres correspondants dans le système de gestion de données de chromatographie, par ex. NeedleChanged et NeedleSeatChanged.

7.10 Entretien de la tête du dispositif de mesure

7.10.1 Composant de la tête du dispositif de mesure



Illustration 158: Vue d'ensemble de la tête du dispositif de mesure

N°	Description
1	Tête du dispositif de mesure
2	Vis de fixation
3	Sortie de rinçage des joints
4	Capillaire vers le capteur de pression
5	Capillaire vers la vanne d'injection
6	Entrée de rinçage des joints

7.10.2 Retrait et raccordement des capillaires des têtes des dispositifs de mesure

- 1. Préparez le passeur d'échantillon pour la maintenance :
 - Arrêtez le débit de la pompe. Patientez jusqu'à ce que la pression du système redescende à zéro avant de poursuivre l'arrêt du passeur d'échantillon.
 - Pour vider la pompe d'évacuation, laissez-la allumée jusqu'à ce qu'il n'y a plus d'eau de condensation dans la conduite d'évacuation. Ensuite, mettez la pompe d'évacuation hors tension et retirez le tube péristaltique de la pompe d'évacuation. Tout liquide restant présent dans le tube de la pompe d'évacuation s'écoule vers l'évacuation.
 - Installez l'unité d'aiguille en position d'entretien et verrouillez-la en place. Reportez-vous à la section Verrouillage de l'unité d'aiguille (> page 251).

Déconnexion des capillaires de la tête du dispositif de mesure

- 1. Déconnectez le capillaire de la vanne d'injection de l'orifice frontal de la tête du dispositif de mesure.
- 2. Déconnectez le capillaire du capteur de pression de l'orifice inférieur de la tête du dispositif de mesure.
- Si vous comptez reconnecter les mêmes capillaires par la suite, installez des capuchons protecteurs sur les extrémités ouvertes des capillaires.

Connexion des capillaires de la tête du dispositif de mesure

- 1. Raccordez le capillaire de la vanne d'injection à l'orifice frontal de la tête du dispositif de mesure.
- Connectez le capillaire du capteur de pression de l'orifice inférieur de la tête du dispositif de mesure.
 Pour un accès plus facile, les vis de fixation des capteurs de pression peuvent être desserrées puis resserrées.
- 3. Finalisez la maintenance du passeur d'échantillon :
 - Déverrouillez l'unité d'aiguille. Reportez-vous à la section Déverrouillage de l'unité d'aiguille (> page 252).
 - Redémarrez le passeur d'échantillon.
 - Préparez le passeur d'échantillon pour l'utilisation. Reportezvous à la section Préparer le système à l'utilisation (> page 161).

7.10.3 Remplacement de la tête du dispositif de mesure

À quel moment exécuter cette opération ?

Tête du dispositif de mesure endommagée

Éléments nécessaires

Tête du dispositif de mesure de remplacement

Outils nécessaires

- Tournevis type Torx T30
- Brucelles (facultatif)

Préparatifs

- 1. Exécutez le script : **C22 Remplacer la tête de mesure** pour déverrouiller la tête du dispositif de mesure.
- 2. Préparez le passeur d'échantillon pour la maintenance :
 - Arrêtez le débit de la pompe. Patientez jusqu'à ce que la pression du système redescende à zéro avant de poursuivre l'arrêt du passeur d'échantillon.
 - Pour vider la pompe d'évacuation, laissez-la allumée jusqu'à ce qu'il n'y a plus d'eau de condensation dans la conduite d'évacuation. Ensuite, mettez la pompe d'évacuation hors tension et retirez le tube péristaltique de la pompe d'évacuation. Tout liquide restant présent dans le tube de la pompe d'évacuation s'écoule vers l'évacuation.
 - Installez l'unité d'aiguille en position d'entretien et verrouillez-la en place. Reportez-vous à la section Verrouillage de l'unité d'aiguille (> page 251).
- 3. Videz le système de rinçage des joints :
 - a) Dévissez le bouchon du réservoir de liquide de rinçage des joints.
 - b) Retirez la conduite de rinçage des joints en même temps que le bouchon du réservoir de liquide de rinçage des joints.
 - c) Pour vider la conduite de rinçage, soulevez le levier de la pompe de rinçage des joints. Relâchez le levier lorsque la conduite est vide.
- Déconnectez les capillaires de la tête du dispositif de mesure. Reportez-vous à la section Retrait et raccordement des capillaires des têtes des dispositifs de mesure (▶ page 263).
- Déconnectez les conduites de rinçage des joints de la tête du dispositif de mesure. Pour faciliter leur retrait, déconnectez les raccords de tube de la tête du dispositif de mesure en laissant les conduites raccordées, et en vous aidant des brucelles.

Retrait de la tête du dispositif de mesure :

- 1. À l'aide du tournevis, desserrez les deux vis qui fixent la tête du dispositif de mesure au passeur d'échantillon.
- 2. Retirez la tête du dispositif de mesure :
 - a) Saisissez la tête du dispositif de mesure par sa face avant. Déplacez délicatement la tête du dispositif de mesure à gauche et à droite plusieurs fois pour détacher le piston du dispositif de mesure.
 - b) Assurez-vous que la tête du dispositif de mesure dépasse d'environ 5 mm du panneau avant du passeur d'échantillon.
 - c) Abaissez avec prudence l'avant de la tête du dispositif de mesure et tirez sur la tête du dispositif de mesure pour le dégager totalement du panneau avant.
 - d) Vérifiez que la tête du dispositif de mesure est retirée avec le piston.



Illustration 159: Retrait de la tête du dispositif de mesure

N°	Description
1	Piston
2	Tête du dispositif de mesure

NOTE Si le piston reste coincé dans le dispositif de mesure, retirez le piston avant de poursuivre la procédure de remplacement. Reportezvous à la section.

Installation de la tête du dispositif de mesure

- Déballez la tête du dispositif de mesure de remplacement. Laissez les bouchons de raccordement installés et les conduites de rinçage des joints raccordées sur la tête du dispositif de mesure de remplacement.
- 2. Introduisez la tête du dispositif de mesure dans l'ouverture du panneau avant.
- 3. Serrez au maximum les deux vis situées sur la tête du dispositif de mesure à l'aide du tournevis (Torx T30).
- 4. Retirez les bouchons de raccordement de la tête du dispositif de mesure.
- 5. Déconnectez les conduites de rinçage des joints d'interconnexion du raccord de tube d'interconnexion. Les conduites de rinçage des joints peuvent rester raccordées à la tête du dispositif de mesure.
- 6. Connectez les conduites de rinçage des joints de la nouvelle tête du dispositif de mesure au système de rinçage des joints. Suivez les étapes d'installation correspondantes pour les conduites de rinçage des joints figurant à la section Remplacement des conduites de rinçage des joints
- Branchez les capillaires sur la tête du dispositif de mesure. Reportezvous à la section Retrait et raccordement des capillaires des têtes des dispositifs de mesure (▶ page 263).
- 8. Finalisez la maintenance du passeur d'échantillon :
 - Déverrouillez l'unité d'aiguille. Reportez-vous à la section Déverrouillage de l'unité d'aiguille (▶ page 252).
 - Redémarrez le passeur d'échantillon.
 - Préparez le passeur d'échantillon pour l'utilisation. Reportezvous à la section Préparer le système à l'utilisation (> page 161).
- Exécutez le script : C22 Remplacer la tête de mesure. Le passeur d'échantillon prépare la tête du dispositif de mesure pour son utilisation.
- Vérifiez s'il y a des fuites dans le système de rinçage des joints. Reportez-vous à la section Test d'étanchéité du système de rinçage des joints () page 282).
- Rincez le système de rinçage des joints. Exécutez le script B01 Modifier les liquides / solvants (avec : Liquide de rinçage des joints arrière et Rafraîchir Uniquement = désactivé.)

7.11 Entretien de la pompe d'évacuation

La pompe d'évacuation (pompe péristaltique) élimine toute condensation qui a pu se créer compartiment à échantillons lors du refroidissement des échantillons et l'évacue via un orifice d'évacuation distinct.



Illustration 160: Pompe d'évacuation

N°	Description
1	Levier de la pompe
2	Rotor de la pompe d'évacuation
3	Tube péristaltique
4	Support de tube
5	Tube de condensation, vers le compartiment à échantillons
6	Tube d'évacuation, vers l'orifice d'évacuation

7.11.1 Test des tubes de la pompe d'évacuation

Testez les tubes de la pompe d'évacuation pour voir s'ils sont obstrués ou s'ils fuient.

À quel moment exécuter cette opération ?

À intervalles réguliers

Éléments nécessaires

Seringue (12 mL)

Préparatifs

- 1. Arrêtez la pompe d'évacuation.
- 2. Aspirez de l'air dans la seringue.

3. Déconnectez le tube d'évacuation et le raccord de tube coudé du tube péristaltique.



Illustration 161: Déconnexion du tube d'évacuation et du raccord de tube coudé

- 4. Introduisez la seringue dans l'extrémité ouverte du tube péristaltique.
- 5. Pour permettre un bon écoulement de l'air dans tous les tubes de la pompe d'évacuation, poussez le levier de la pompe d'évacuation vers le haut, puis maintenez-le d'une main dans cette position.



Illustration 162: Levier de la pompe d'évacuation poussé vers le haut

Recherche d'obstruction

- 1. Appuyez sur le piston de la seringue pour injecter l'air dans le tube.
- 2. Déterminez si vous ressentez une résistance lorsque vous enfoncez le piston de la seringue :

Observation	Étapes
ll y a une résistance	Le tube péristaltique peut être obstrué. Retirez la seringue et remplacez le tube péristaltique (reportez-vous à la section Remplacement du tube de la pompe d'évacuation () page 270)).
Il n'y a aucune résistance	Recherchez des fuites.

Recherche de fuites

1. Aspirez de l'air dans la seringue.

2. Déconnectez le tube de condensation et le raccord de tube coudé du tube péristaltique.



Illustration 163: Déconnexion du tube de condensation et du raccord de tube coudé

- 3. Obstruez une extrémité du tube péristaltique à l'aide d'un doigt.
- 4. Appuyez sur le piston de la seringue pour injecter l'air dans le tube.
- 5. Déterminez si vous ressentez que la pression augmente lorsque vous enfoncez le piston de la seringue :

Observation	Étapes
Il n'y a aucune pression	Le tube péristaltique peut avoir une fuite ou être endommagé. Retirez la seringue et remplacez le tube péristaltique (reportez-vous à la section Remplacement du tube de la pompe d'évacuation () page 270)).
Il y a de la pression	Retirez la seringue et reconnectez les raccords de tube coudés, le tube de condensation et le tube d'évacuation. Relâchez le levier de la pompe.

7.11.2 Remplacement du tube de la pompe d'évacuation

À quel moment exécuter cette opération ?

- Tous les six mois en tant que procédure de maintenance préventive pour la pompe d'évacuation
- Le tube de la pompe d'évacuation est obstrué (reportez-vous à la section Test des tubes de la pompe d'évacuation (> page 268))
- Le tube de la pompe d'évacuation a une fuite ou est endommagé (reportez-vous à la section Test des tubes de la pompe d'évacuation (> page 268))

Éléments nécessaires

- Tube péristaltique de remplacement pour la pompe d'évacuation
- 2 raccords de tube coudés pour le tube de condensation et le tube d'évacuation

Outils nécessaires

Coupe-tube (en option)

Préparatifs

- 1. Arrêtez la pompe d'évacuation.
- Déconnectez le tube d'évacuation et le tube de condensation des raccords de tube sur le tube de la pompe d'évacuation (tube péristaltique). Laissez les raccords de tube connectés au tube péristaltique.



Illustration 164: Déconnexion du tube d'évacuation et du tube de condensation

- 3. Déballez le tube péristaltique de remplacement et les raccords de tube coudés.
- 4. Connectez les raccords de tube coudés à chaque extrémité du tube péristaltique de remplacement.

Procédez comme suit

1. Actionnez vers le haut le levier de la pompe d'évacuation, puis maintenez-le d'une main dans cette position.



Illustration 165: Levier de pompe poussé vers le haut

- 2. Avec l'autre main, dégagez le tube péristaltique du rotor de la pompe d'évacuation, puis retirez-le de la pompe d'évacuation. Relâchez le levier de la pompe.
- Vérifiez que le tube péristaltique de remplacement est de la même longueur que celui installé précédemment. Si nécessaire, ajustez la longueur du tube péristaltique de remplacement à l'aide du coupetube.
- Enfoncez le tube péristaltique de remplacement dans les supports de tube. Vérifiez que le tube est correctement inséré dans les supports.
- Actionnez vers le haut le levier de la pompe d'évacuation. Insérez le tube péristaltique de rechange entre le levier de la pompe et le rotor de la pompe d'évacuation, puis enroulez-le autour du rotor.
- 6. Relâchez le levier de la pompe.
- 7. Connectez le tube d'évacuation et le tube de condensation aux raccords de tube sur le tube péristaltique.



Illustration 166: Connexion du tube d'évacuation et du tube de condensation

8. Mettez la pompe d'évacuation sous tension pour vérifier son bon fonctionnement.

7.12 Entretien du dispositif de rinçage de l'aiguille

Pour un aperçu de l'ensemble des conduites de rinçage, reportez-vous à la section Dispositif de rinçage de l'aiguille (vue d'ensemble) (▶ page 58).

7.12.1 Remplacement des conduites de rinçage entre le réservoir et l'orifice d'admission de la pompe de rinçage de l'aiguille

Suivez les étapes ci-dessous pour remplacer les deux conduites de rinçage entre les réservoirs de rinçage de l'aiguille et les orifices d'admission de la pompe de rinçage de l'aiguille du passeur d'échantillon.

À quel moment exécuter cette opération ?

La conduite de rinçage est obstruée ou endommagée

Pièces et éléments complémentaires nécessaires

- Liquide de rinçage de l'aiguille fraîchement préparé
- Remplacement des conduites de rinçage étiquetées SWP (rouge) et WWP (verte)

Outils nécessaires

Coupe-tube (en option)

NOTE 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 de la conduite.

Préparatifs

- 1. Préparez du liquide de rinçage de l'aiguille frais.
- 2. Videz le dispositif de rinçage de l'extérieur de l'aiguille :
 - a) Dévissez les bouchons des réservoirs de rinçage de l'aiguille.
 - b) Retirez les conduites de rinçage (étiquetées WWP et SWP) du réservoir en même temps que le bouchon du réservoir.
 - c) Lorsque vous retirez les conduites des guides de tubulure, veillez à ne pas enlever d'autres tubulures présentes dans les guides.
 - d) Placez les conduites de rinçage dans un récipient pour déchets pour les vider.

Procédez comme suit

- 1. Retirez le guide de maintien des conduites de rinçage.
- 2. Sortez les conduites des bouchons de réservoir.
- 3. Déconnectez les conduites de rinçage des orifices d'aspiration de la pompe de rinçage de l'aiguille.



Illustration 167: Déconnexion des conduites de rinçage de l'orifice d'admission de la pompe de rinçage de l'aiguille

- 4. Connectez les conduites de rinçage de rechange aux orifices d'admission de la pompe de rinçage de l'aiguille :
 - a) Enfoncez le tube rigide dans le connecteur.
 - b) Poussez le tube en silicone par dessus le connecteur.
- Acheminez les conduites de rinçage depuis le passeur d'échantillon jusqu'au bac à solvants, en passant par les guides de tubulure dans le passeur d'échantillon et dans tous les modules situés au-dessus du passeur d'échantillon dans le système empilé.
- Acheminez les conduites de rinçage à travers le trou de guidage situé à l'intérieur du bac à solvants et poussez-les dans un ou plusieurs guides de tubulure.



Illustration 168: Trou de guidage et guides de tubulure dans le bac à solvants

N°	Description
А	Trou de guidage (bac à solvants)
1-4	Guides de tubulure

- 7. Faites passer chaque conduite de rinçage de l'aiguille par un guide de maintien.
- 8. Faites passer chaque conduite de rinçage de l'aiguille par un trou du bouchon du réservoir de rinçage. Le guide de maintien permet d'immobiliser la conduite de solvant dans le réservoir.
- 9. Obturez les trous restants des bouchons de réservoir au moyen de capuchons protecteurs.
- 10. Glissez un porte-filtre avec adaptateur sur chaque conduite de rinçage.
- Remplissez les réservoirs de rinçage de l'aiguille avec du liquide de rinçage de l'aiguille fraîchement préparé et redémarrez le dispositif de rinçage de l'aiguille. Pour en savoir plus, consultez la section Remplissage des réservoirs de rinçage de l'aiguille (▶ page 131).

7.12.2 Remplacement de la conduite de rinçage entre l'orifice de refoulement de la pompe de rinçage de l'aiguille et l'orifice de rinçage

Suivez les étapes ci-dessous pour remplacer la conduite de rinçage entre les orifices de refoulement de la pompe de rinçage de l'aiguille et l'orifice de rinçage de l'aiguille.

À quel moment exécuter cette opération ?

La conduite de rinçage est obstruée ou endommagée

Pièces et éléments complémentaires nécessaires

- Liquide de rinçage de l'aiguille fraîchement préparé
- Remplacement de la conduite de rinçage

Préparatifs

- 1. Préparez du liquide de rinçage de l'aiguille frais.
- 2. Videz le dispositif de rinçage de l'extérieur de l'aiguille :
 - a) Dévissez les bouchons des réservoirs de rinçage de l'aiguille.
 - b) Retirez les conduites de rinçage (étiquetées WWP et SWP) du réservoir en même temps que le bouchon du réservoir.
 - c) Lorsque vous retirez les conduites des guides de tubulure, veillez à ne pas enlever d'autres tubulures présentes dans les guides.
 - d) Placez les conduites de rinçage dans un récipient pour déchets pour les vider.

Procédez comme suit

1. Déconnectez les deux extrémités de la conduite de rinçage des orifices de refoulement de la pompe de rinçage de l'aiguille.



Illustration 169: Déconnexion des conduites de rinçage de l'orifice de refoulement de la pompe de rinçage de l'aiguille

2. Déconnectez la conduite de rinçage de l'orifice d'admission de rinçage.



Illustration 170: Déconnexion de la conduite de rinçage de l'orifice de rinçage

- Connectez la conduite de rinçage de rechange aux orifices de refoulement de la pompe de rinçage de l'aiguille et à l'orifice d'admission de rinçage.
- Remplissez les réservoirs de rinçage de l'aiguille avec du liquide de rinçage de l'aiguille fraîchement préparé et redémarrez le dispositif de rinçage de l'aiguille. Pour en savoir plus, consultez la section Remplissage des réservoirs de rinçage de l'aiguille (▶ page 131).

7.12.3 Remplacement des conduites de rinçage entre les orifices d'admission du réservoir et du dégazeur

Suivez les étapes ci-dessous pour remplacer les deux conduites de rinçage entre les réservoirs de rinçage de l'aiguille et l'orifice d'admission du dégazeur placé dans la pompe.

À quel moment exécuter cette opération ?

La conduite de rinçage entre le réservoir et le dégazeur est obstruée ou endommagée

Pièces et éléments complémentaires nécessaires

- Liquide de rinçage de l'aiguille fraîchement préparé
- Remplacement des conduites de rinçage étiquetées W (verte) et S (rouge)

Outils nécessaires

Coupe-tube (en option)

NOTE 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 de la conduite.

Préparatifs

- 1. Préparez du liquide de rinçage de l'aiguille frais.
- 2. Videz le dispositif de rinçage de l'aiguille :
 - a) Dévissez les bouchons des réservoirs de rinçage de l'aiguille.
 - b) Retirez les conduites de rinçages (étiquetées W et S) des réservoirs en même temps que le bouchon.
 - c) Placez les conduites de rinçage dans un récipient pour déchets pour les vider.

Procédez comme suit

- 1. Retirez le guide de maintien des conduites de rinçage.
- 2. Sortez les conduites des bouchons de réservoir.
- 3. Déconnectez les conduites de rinçage du dégazeur.

4. Connectez les conduites de rinçage de rechange aux orifices d'admission du dégazeur.





N°	Description
1	Orifice d'admission pour la conduite de rinçage étiquetée W (verte) vers le réservoir
2	Orifice d'admission pour la conduite de rinçage étiquetée S (rouge) vers le réservoir
3	Orifice de refoulement pour la conduite de rinçage étiquetée S (rouge) vers le passeur d'échantillon
4	Orifice de refoulement pour la conduite de rinçage étiquetée W (verte) vers le passeur d'échantillon

5. Acheminez les conduites de rinçage depuis le passeur d'échantillon jusqu'au bac à solvants, en passant par les guides de tubulure dans le passeur d'échantillon et dans tous les modules situés au-dessus du passeur d'échantillon dans le système empilé.

6. Acheminez les conduites de rinçage à travers le trou de guidage situé à l'intérieur du bac à solvants et poussez-les dans un ou plusieurs guides de tubulure.



Illustration 172: Trou de guidage et guides de tubulure dans le bac à solvants

N°	Description
А	Trou de guidage (bac à solvants)
1-4	Guides de tubulure

- 7. Faites passer chaque conduite de rinçage de l'aiguille par un guide de maintien.
- Faites passer chaque conduite de rinçage de l'aiguille par un trou du bouchon du réservoir de rinçage. Le guide de maintien permet d'immobiliser la conduite de solvant dans le réservoir.
- 9. Obturez les trous restants des bouchons de réservoir au moyen de capuchons protecteurs.
- 10. Glissez un porte-filtre avec adaptateur sur chaque conduite de rinçage.
- 11. Remplissez les réservoirs de rinçage de l'aiguille avec du liquide de rinçage de l'aiguille fraîchement préparé et redémarrez le dispositif de rinçage de l'aiguille. Reportez-vous à la section Remplissage des réservoirs de rinçage de l'aiguille (▶ page 131).

7.12.4 Remplacement des conduites de rinçage entre l'orifice de refoulement du dégazeur vers la vanne de sélection

Suivez les étapes ci-dessous pour remplacer les deux conduites de rinçage entre l'orifice de refoulement du dégazeur placé dans la pompe et les orifices d'admission de rinçage de l'aiguille de la vanne de sélection du passeur d'échantillon.

À quel moment exécuter cette opération ?

La conduite de rinçage est obstruée ou endommagée

Pièces et éléments complémentaires nécessaires

- Liquide de rinçage de l'aiguille fraîchement préparé
- Remplacement des conduites de rinçage

Préparatifs

- 1. Préparez du liquide de rinçage de l'aiguille frais.
- 2. Videz le dispositif de rinçage de l'aiguille :
 - a) Dévissez les bouchons des réservoirs de rinçage de l'aiguille.
 - b) Retirez les conduites de rinçages (étiquetées W et S) des réservoirs en même temps que le bouchon.
 - c) Placez les conduites de rinçage dans un récipient pour déchets pour les vider.

Procédez comme suit

- 1. Retirez le guide de maintien des conduites de rinçage.
- 2. Déconnectez les conduites de rinçage de la vanne de sélection du passeur d'échantillon.
- 3. Déconnectez les conduites de l'orifice de refoulement du dégazeur.
- 4. Connectez les conduites de rinçage de rechange aux orifices de refoulement du dégazeur.
- 5. Connectez les conduites de rinçage de l'aiguille à la vanne de sélection du passeur d'échantillon.
- 6. Remplissez les réservoirs de rinçage de l'aiguille avec du liquide de rinçage de l'aiguille fraîchement préparé et redémarrez le dispositif de rinçage de l'aiguille. Reportez-vous à la section Remplissage des réservoirs de rinçage de l'aiguille (▶ page 131).

Consultez également

Dispositif de rinçage de l'aiguille (vue d'ensemble) (> page 58)



7.12.5 Remplacement du clapet de non-retour du rinçage de l'aiguille

Illustration 173: Vanne de sélection avec clapet de non-retour

N°	Description
1	Capillaire Viper vers la vanne de solvant
2	Clapet de non-retour pour rinçage de l'aiguille
3	Vanne de sélection pour rinçage de l'aiguille

À quel moment exécuter cette opération ?

Le clapet de non-retour est défectueux

Éléments nécessaires

Clapet de non-retour de remplacement

Procédez comme suit

- 1. Desserrez le connecteur Viper sur le clapet de non-retour et déconnectez le capillaire Viper de la vanne de solvant.
- 2. Dévissez le clapet de non-retour de la vanne de sélection.
- 3. Installez le clapet de non-retour de rechange.
- 4. Connectez le capillaire Viper à la vanne de solvant. Serrez le raccord.
- 5. Purgez le passeur d'échantillon.

Consultez également

- Attributions des orifices de la vanne du passeur d'échantillon (> page 48)
- Purge (vue d'ensemble) (▶ page 65)

7.13 Entretien du système de rinçage des joints

La présente section décrit comment entretenir et remplacer les éléments du système de rinçage des joints. Suivez les instructions de la procédure de maintenance que vous souhaitez effectuer.

7.13.1 Test d'étanchéité du système de rinçage des joints

À quel moment exécuter cette opération ?

- À intervalles réguliers
- Quand vous avez installé la tête de pompe pendant les procédures de maintenance
- Après avoir remplacé la tête du dispositif de mesure

Matériel nécessaire

- Seringue (12 mL)
- le liquide de rinçage des joints ;

Procédez comme suit

1. Débranchez le tube du détecteur de rinçage des joints.





N°	Description
1	Détecteur de rinçage des joints
2	Orifice de raccordement de tube (orifice étiqueté in)

2. Introduisez une seringue dans l'extrémité ouverte du tube.

- Pour permettre un bon écoulement du liquide dans l'ensemble du système de rinçage des joints, soulevez le levier de la pompe de rinçage des joints, puis maintenez-le d'une main dans cette position.
- 4. Tirez sur le piston de la seringue pour aspirer le liquide de rinçage des joints dans la seringue.
- 5. Lorsque la seringue contient environ 10 mL de liquide, relâchez le levier de la pompe de rinçage des joints.
- Levez la conduite du réservoir de liquide de rinçage des joints audessus de la tête du dispositif de mesure et pincez le tube pour le bloquer.
- 7. Appuyez légèrement sur le piston de la seringue pour appliquer de la pression sur le système de rinçage des joints.



Illustration 175: Position dans laquelle tenir la conduite du réservoir de liquide de rinçage des joints

 Vérifiez si du liquide s'échappe des raccordements du système de rinçage des joints, sous les têtes de pompe ou sous la tête du dispositif de mesure.

Observation	Mesure corrective
Un raccordement fuit	 Resserrez ou remplacez les raccordements qui fuient. Répétez le test.
ll y a une fuite sous les têtes de pompe	 Vérifiez si les joints de la tête de pompe et le joint de rinçage des joints sont correctement installés. Répétez le test.
Une fuite est présente sous la tête du dispositif de mesure	 Vérifiez que la tête du dispositif de mesure est bien installée. S'il y a toujours une fuite, remplacez la tête du dispositif de mesure. Répétez le test.
Aucune fuite n'est observée	Retirez la seringue et reconnectez le tube sur le détecteur de rinçage des joints (orifice étiqueté in).

Consultez également

Remplacement de la tête du dispositif de mesure (> page 264)

7.13.2 Remplacement de la conduite de rinçage du réservoir de liquide de rinçage des joints

Suivez les étapes ci-dessous pour remplacer la conduite de rinçage entre le réservoir de liquide de rinçage des joints et l'orifice d'admission de rinçage des joints sur le passeur d'échantillon.



Illustration 176: Position de la conduite du réservoir de liquide de rinçage des joints

À quel moment exécuter cette opération ?

La conduite du réservoir de liquide de rinçage des joints est obstruée ou endommagée

Éléments nécessaires

Tubes en silicone et raccords

Outils nécessaires

Coupe-tube

Préparatifs

- 1. Videz le système de rinçage des joints :
 - a) Dévissez le bouchon du réservoir de liquide de rinçage des joints.
 - b) Retirez la conduite de rinçage des joints en même temps que le bouchon du réservoir de liquide de rinçage des joints.
 - c) Pour vider la conduite de rinçage, soulevez le levier de la pompe de rinçage des joints. Relâchez le levier lorsque la conduite est vide.

Retrait de la conduite du réservoir de liquide de rinçage des joints

1. Tirez sur la conduite du réservoir de liquide de rinçage des joints munie du raccord de tube pour la sortir de l'orifice d'admission de rinçage des joints.



Illustration 177: Sortie de la conduite de rinçage des joints munie du raccord de tube de l'orifice d'admission de rinçage des joints

 Déconnectez la conduite du réservoir de liquide de rinçage des joints, le raccord de tube et la conduite d'admission de rinçage des joints. La conduite de rinçage des joints sur la tête du dispositif de mesure peut rester connectée.



Illustration 178: Déconnexion de la conduite du réservoir de liquide de rinçage des joints et de la conduite d'admission de rinçage des joints

- 3. Déconnectez la conduite du réservoir de liquide de rinçage des joints :
 - a) Retirez le guide de maintien de la conduite de rinçage.
 - b) Sortez la conduite de rinçage du bouchon du réservoir.



Illustration 179: Conduite de rinçage et bouchon du réservoir

N°	Description
1	Conduite de rinçage
2	Guide de maintien
3	Bouchon de réservoir
4	Capuchons protecteurs

Connexion de la nouvelle conduite du réservoir de liquide de rinçage des joints

- 1. Coupez le tube à la longueur requise, le cas échéant.
- Connectez la conduite de rinçage des joints à la conduite d'admission de rinçage des joints à l'aide d'un nouveau raccord de tube.
- 3. Enfoncez le raccord de tube avec les conduites attachées dans l'orifice d'admission de rinçage des joints.
- 4. Acheminez la conduite de rinçage des joints depuis le passeur d'échantillon jusqu'au bac à solvants, en passant par les guides de tubulure dans le passeur d'échantillon et dans tous les modules situés au-dessus du passeur d'échantillon dans le système empilé.
- 5. Acheminez la conduite de rinçage des joints à travers le trou de guidage situé à l'intérieur du bac à solvants et fixez-la dans l'un des guides de tubulure.



Illustration 180: Trou de guidage et guides de tubulure dans le bac à solvants

N°	Description
А	Trou de guidage (bac à solvants)
1-4	Guides de tubulure

6. Raccordez le réservoir de rinçage des joints à la conduite :

- a) Rincez soigneusement le réservoir de rinçage des joints avec un solvant ultrapur.
- b) Remplissez le réservoir de rinçage des joints avec du liquide de rinçage des joints fraîchement préparé.
- c) Faites passer la conduite de rinçage des joints par le guide de maintien et un trou du bouchon du réservoir de rinçage des joints. Le guide de maintien permet d'immobiliser la tubulure dans le réservoir. Obturez les trous restants du bouchon de réservoir au moyen de capuchons protecteurs.

Redémarrage du système de rinçage des joints

- Resserrez le bouchon du réservoir manuellement. Le guide de maintien doit rester dans le trou du bouchon de réservoir. Si ce n'est pas le cas, enfoncez le guide de maintien dans le trou afin d'immobiliser la tubulure dans le bouchon.
- 2. Inspectez les conduites de rinçage sur l'ensemble du trajet d'écoulement :
 - Assurez-vous que les conduites de rinçage ne sont, à aucun endroit du trajet d'écoulement, pliées (plicaturées), pincées ou serrées.
 - b) Si vous devez raccourcir la longueur du tube, utilisez un coupetube. Veillez à ce que la coupe soit à angle droit par rapport à l'axe de la conduite.
- Rincez le système de rinçage des joints. Exécutez le script B01 Modifier les liquides / solvants (avec : Liquide de rinçage des joints arrière et Rafraîchir Uniquement = désactivé.)

Consultez également

Système de rinçage des joints (vue d'ensemble) (page 60)

7.13.3 Remplacement de la conduite d'admission de rinçage des joints

Suivez les étapes ci-dessous pour remplacer la conduite de rinçage entre l'orifice d'admission de rinçage des joints sur le passeur d'échantillon et l'orifice d'admission de rinçage des joints de la tête du dispositif de mesure.



Illustration 181: Position de la conduite d'admission de rinçage des joints

À quel moment exécuter cette opération ?

La conduite d'admission de rinçage des joints est obstruée ou endommagée

Éléments nécessaires

Tubes en silicone et raccords

Outils nécessaires

Coupe-tube

Préparatifs

- 1. Videz le système de rinçage des joints :
 - a) Dévissez le bouchon du réservoir de liquide de rinçage des joints.
 - b) Retirez la conduite de rinçage des joints en même temps que le bouchon du réservoir de liquide de rinçage des joints.
 - c) Pour vider la conduite de rinçage, soulevez le levier de la pompe de rinçage des joints. Relâchez le levier lorsque la conduite est vide.
Procédez comme suit

 Tirez sur la conduite d'admission de rinçage des joints munie du raccord de tube pour la sortir de l'entrée de rinçage des joints sur le haut de la tête du dispositif de mesure.





2. Tirez sur la conduite de rinçage des joints munie du raccord de tube pour la sortir de l'orifice d'admission de rinçage des joints.



Illustration 183: Sortie de la conduite de rinçage des joints munie du raccord de tube de l'orifice d'admission de rinçage des joints

3. Déconnectez la conduite d'admission de rinçage des joints du raccord de tube. La conduite du réservoir de liquide de rinçage des joints, de l'autre côté du raccord de tube, peut rester connectée.



Illustration 184: Déconnexion de la conduite d'admission de rinçage des joints de l'orifice d'admission de rinçage des joints

- 4. Connectez la nouvelle conduite de rinçage des joints à la tête du dispositif de mesure à l'aide d'un nouveau raccord de tube. Coupez le tube à la longueur requise, le cas échéant.
- 5. Dirigez la conduite d'admission de rinçage des joints jusqu'à l'orifice d'admission de rinçage des joints.
- 6. Connectez la conduite d'admission de rinçage des joints au raccord de tube.
- 7. Enfoncez le raccord de tube avec les conduites attachées dans l'orifice d'admission de rinçage des joints.

Redémarrage du système de rinçage des joints

- Resserrez le bouchon du réservoir manuellement. Le guide de maintien doit rester dans le trou du bouchon de réservoir. Si ce n'est pas le cas, enfoncez le guide de maintien dans le trou afin d'immobiliser la tubulure dans le bouchon.
- 2. Inspectez les conduites de rinçage sur l'ensemble du trajet d'écoulement :
 - Assurez-vous que les conduites de rinçage ne sont, à aucun endroit du trajet d'écoulement, pliées (plicaturées), pincées ou serrées.
 - b) Si vous devez raccourcir la longueur du tube, utilisez un coupetube. Veillez à ce que la coupe soit à angle droit par rapport à l'axe de la conduite.
- Rincez le système de rinçage des joints. Exécutez le script B01 Modifier les liquides / solvants (avec : Liquide de rinçage des joints arrière et Rafraîchir Uniquement = désactivé.)

7.13.4 Remplacement de la conduite d'admission de rinçage des joints de la pompe

Suivez les étapes ci-dessous pour remplacer la conduite de rinçage entre l'orifice de refoulement de la tête du dispositif de mesure et l'orifice d'admission de la pompe de rinçage des joints.



Illustration 185: Position de la conduite d'admission de la pompe de rinçage des joints

À quel moment exécuter cette opération ?

La conduite d'admission de la pompe de rinçage des joints est obstruée ou endommagée

Éléments nécessaires

Tubes en silicone et raccords

Outils nécessaires

Coupe-tube

Préparatifs

- 1. Videz le système de rinçage des joints :
 - a) Dévissez le bouchon du réservoir de liquide de rinçage des joints.
 - b) Retirez la conduite de rinçage des joints en même temps que le bouchon du réservoir de liquide de rinçage des joints.
 - c) Pour vider la conduite de rinçage, soulevez le levier de la pompe de rinçage des joints. Relâchez le levier lorsque la conduite est vide.

Procédez comme suit

1. Tirez sur le raccord de tube avec la conduite de la pompe de rinçage des joints pour le sortir de l'orifice de refoulement de rinçage des joints sur la tête du dispositif de mesure.



Illustration 186: Déconnexion de la conduite d'admission de la pompe de rinçage des joints de la tête du dispositif de mesure

2. Déconnectez la conduite d'admission de la pompe de rinçage des joints du tube péristaltique sur la pompe de rinçage des joints dans la pompe.



Illustration 187: Déconnexion de la conduite d'admission de la pompe de rinçage des joints

- 3. Connectez un nouveau raccord de tube sur la nouvelle conduite d'admission de la pompe de rinçage des joints. Coupez le tube à la longueur requise, le cas échéant.
- 4. Connectez la nouvelle conduite d'admission de la pompe de rinçage des joints à la tête du dispositif de mesure.
- 5. Connectez la conduite d'admission de la pompe de rinçage des joints au tube péristaltique de la pompe de rinçage des joints.

Redémarrage du système de rinçage des joints

- Resserrez le bouchon du réservoir manuellement. Le guide de maintien doit rester dans le trou du bouchon de réservoir. Si ce n'est pas le cas, enfoncez le guide de maintien dans le trou afin d'immobiliser la tubulure dans le bouchon.
- 2. Inspectez les conduites de rinçage sur l'ensemble du trajet d'écoulement :
 - Assurez-vous que les conduites de rinçage ne sont, à aucun endroit du trajet d'écoulement, pliées (plicaturées), pincées ou serrées.
 - b) Si vous devez raccourcir la longueur du tube, utilisez un coupetube. Veillez à ce que la coupe soit à angle droit par rapport à l'axe de la conduite.
- Rincez le système de rinçage des joints. Exécutez le script B01 Modifier les liquides / solvants (avec : Liquide de rinçage des joints arrière et Rafraîchir Uniquement = désactivé.)

7.13.5 Remplacement du tube péristaltique dans la pompe de rinçage des joints

Suivez les étapes ci-dessous pour remplacer le tube péristaltique dans la pompe de rinçage des joints.



Illustration 188: Position du tube péristaltique dans la pompe de rinçage des joints

À quel moment exécuter cette opération ?

- Le tube est obstrué ou endommagé
- Tous les 6 mois en tant que procédure de maintenance préventive pour la pompe de rinçage des joints

Éléments nécessaires

Tube péristaltique (tube PharMed[™]), raccords de tube et supports

Outils nécessaires

Coupe-tube

Préparatifs

- 1. Videz le système de rinçage des joints :
 - a) Dévissez le bouchon du réservoir de liquide de rinçage des joints.
 - b) Retirez la conduite de rinçage des joints en même temps que le bouchon du réservoir de liquide de rinçage des joints.
 - c) Pour vider la conduite de rinçage, soulevez le levier de la pompe de rinçage des joints. Relâchez le levier lorsque la conduite est vide.

Remplacement du tube péristaltique



Illustration 189: Pompe de rinçage des joints

N°	Description
1	Levier de la pompe (soulevé)
2	Rotor de la pompe
3	Conduite de rinçage des joints du passeur d'échantillon (tubes en silicone)
4	Raccords de tube
5	Supports de tube pour le tube péristaltique
6	Conduite de rinçage des joints vers la tête de pompe (tubes en silicone)
7	Tubes péristaltiques

- 1. Déconnectez le tube péristaltique des raccords de tube qui le connectent au tube en silicone.
- 2. Tirez le levier situé sur la pompe de rinçage des joints vers le haut, et retirez le tube péristaltique de la pompe et des supports de tube.
- Connectez le nouveau tube péristaltique au tube en silicone à l'aide de raccords de tube. Coupez le tube à la longueur requise, le cas échéant.
- 4. Insérez le tube péristaltique dans le support de tube supérieur droit.
- 5. Tirez le levier situé sur la pompe de rinçage des joints vers le haut, puis maintenez-le d'une main dans cette position.
- 6. Avec l'autre main, insérez le tube entre le levier et le rotor, puis enroulez-le autour du rotor.

- 7. Relâchez le levier.
- 8. Insérez le tube péristaltique dans le support de tube inférieur droit.
- 9. Insérez le tube péristaltique dans le support de tube gauche.
- Vérifiez que le tube péristaltique est correctement inséré dans les supports. Dans le cas contraire, enfoncez le tube dans les supports. Assurez-vous que le tube n'est pas pincé ou serré dans les supports.
- 11. Redémarrez le système de rinçage des joints.

Redémarrage du système de rinçage des joints

- Resserrez le bouchon du réservoir manuellement. Le guide de maintien doit rester dans le trou du bouchon de réservoir. Si ce n'est pas le cas, enfoncez le guide de maintien dans le trou afin d'immobiliser la tubulure dans le bouchon.
- 2. Inspectez les conduites de rinçage sur l'ensemble du trajet d'écoulement :
 - Assurez-vous que les conduites de rinçage ne sont, à aucun endroit du trajet d'écoulement, pliées (plicaturées), pincées ou serrées.
 - b) Si vous devez raccourcir la longueur du tube, utilisez un coupetube. Veillez à ce que la coupe soit à angle droit par rapport à l'axe de la conduite.
- Rincez le système de rinçage des joints. Exécutez le script B01 Modifier les liquides / solvants (avec : Liquide de rinçage des joints arrière et Rafraîchir Uniquement = désactivé.)

7.13.6 Remplacement des conduites de rinçage des joints entre la pompe de rinçage des joints et le détecteur de rinçage des joints

Suivez les étapes ci-dessous pour remplacer les conduites de rinçage entre le tube péristaltique (pompe de rinçage des joints) et le détecteur de rinçage des joints.



Illustration 190: Conduites de rinçage des joints entre le tube péristaltique et le détecteur de rinçage des joints

N°	Description
1	Conduite de rinçage entre le tube péristaltique et la tête de pompe gauche
2	Conduite de rinçages sur la tête de pompe gauche
3	Conduite de rinçage entre les têtes de pompe
4	Conduite de rinçage sur la tête de pompe droite
5	Conduite de rinçage entre la tête de pompe droite et le détecteur de rinçage des joints

À quel moment exécuter cette opération ?

Les conduites de rinçage des joints sont obstruées ou endommagées

Éléments nécessaires

Tubulure en silicone

Outils nécessaires

Coupe-tube

Préparatifs

- 1. Videz le système de rinçage des joints :
 - a) Dévissez le bouchon du réservoir de liquide de rinçage des joints.
 - b) Retirez la conduite de rinçage des joints en même temps que le bouchon du réservoir de liquide de rinçage des joints.
 - c) Pour vider la conduite de rinçage, soulevez le levier de la pompe de rinçage des joints. Relâchez le levier lorsque la conduite est vide.

Procédez comme suit

- 1. Déconnectez le tube de rinçage des joints des raccords de tube. Ne retirez pas les raccords de tube.
- 2. Connectez les nouvelles conduites de rinçage des joints. Coupez le tube à la longueur requise, le cas échéant.

Redémarrage du système de rinçage des joints

- Resserrez le bouchon du réservoir manuellement. Le guide de maintien doit rester dans le trou du bouchon de réservoir. Si ce n'est pas le cas, enfoncez le guide de maintien dans le trou afin d'immobiliser la tubulure dans le bouchon.
- 2. Inspectez les conduites de rinçage sur l'ensemble du trajet d'écoulement :
 - Assurez-vous que les conduites de rinçage ne sont, à aucun endroit du trajet d'écoulement, pliées (plicaturées), pincées ou serrées.
 - b) Si vous devez raccourcir la longueur du tube, utilisez un coupetube. Veillez à ce que la coupe soit à angle droit par rapport à l'axe de la conduite.
- Rincez le système de rinçage des joints. Exécutez le script B01 Modifier les liquides / solvants (avec : Liquide de rinçage des joints arrière et Rafraîchir Uniquement = désactivé.)

7.13.7 Remplacement du détecteur de rinçage des joints

À quel moment exécuter cette opération ?

- Contamination des électrodes du détecteur
- Dysfonctionnement du système de rinçage des joints

Éléments nécessaires

Bloc de détecteur de rinçage des joints, incluant le détecteur et le pied

Préparatifs

Pour préparer le nouveau bloc de détecteur de rinçage des joints, enfoncez le détecteur de rinçage des joints dans le pied. Contrôlez l'orientation du pied.



Illustration 191: Montage des pièces du détecteur de rinçage des joints

N°	Description
1a	Pied (bas)
1b	Pied (haut)
2	Détecteur de rinçage des joints

Procédez comme suit



Illustration 192: Retrait du bloc de détecteur de rinçage des joints

N°	Description
1	Pied, détecteur de rinçage des joints
2	Détecteur de rinçage des joints
3	Câble du détecteur
4	Conduite de rinçage des joints

- 1. Déconnectez la conduite de rinçage des joints du détecteur de rinçage des joints (orifice étiqueté **in**).
- 2. Déconnectez le câble du détecteur de l'orifice **DROP DET**.
- 3. Saisissez et tirez le bloc du détecteur de rinçage des joints vers le haut par le pied.
- 4. Poussez le nouveau bloc de détecteur de rinçage des joints dans la pompe.



Illustration 193: Installation du bloc de détecteur de rinçage des joints

- 5. Rebranchez la conduite de rinçage des joints sur le détecteur de rinçage des joints (orifice étiqueté **in**).
- 6. Connectez le câble du détecteur à l'orifice DROP DET.

 Rincez le système de rinçage des joints. Exécutez le script B01 – Modifier les liquides / solvants (avec : Liquide de rinçage des joints arrière et Rafraîchir Uniquement = désactivé.)

7.14 Inspection des vannes du passeur d'échantillon à la recherche de fuites

À quel moment exécuter cette opération ?

- Des gouttelettes de liquide apparaissent à un raccord Viper
- Des gouttelettes de liquide apparaissent au trou d'évacuation



Illustration 194: Trous d'évacuation sur les vannes du passeur d'échantillon

Outils nécessaires

Tournevis dynamométrique avec embout Viper et embout pour le siège de l'aiguille

Procédez comme suit

- Des gouttelettes de liquide apparaissent au trou d'évacuation : Essuyez les gouttelettes et observez la vanne. Si des gouttelettes de liquide continuent à apparaître au trou d'évacuation, prenez contact avec l'assistance technique de Thermo Fisher Scientific. Sinon, poursuivez comme suit.
- 2. Comme les fuites surviennent généralement aux raccords, inspectez visuellement tous les raccords du trajet d'écoulement.
- Desserrez les raccords des capillaires à la vanne touchée et resserrez-les avec le tournevis dynamométrique muni de l'embout Viper.

AVIS

- Serrez ou desserrez les capillaires Viper uniquement avec vos doigts et avec l'outil dynamométrique spécifique. N'utilisez pas d'outils autres que ceux fournis par Thermo Fisher Scientific à cette fin. Utilisez cet outil uniquement pour les raccordements pour lesquels il est prévu.
- 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.
- 4. Si la vanne touchée est la vanne d'injection avec le siège de l'aiguille :
 Desserrez et resserrez le siège de l'aiguille avec le tournevis dynamométrique et l'embout pour le siège de l'aiguille.

AVIS—Le siège de l'aiguille pourrait être endommagé par l'utilisation d'autres outils

Le tournevis dynamométrique avec l'embout pour le siège de l'aiguille est destiné au serrage du siège de l'aiguille. À l'aide de cet outil, le siège de l'aiguille est correctement serré sans serrage excessif. N'utilisez pas un autre outil.

AVIS—Le tournevis dynamométrique avec l'embout pour le siège de l'aiguille peut être endommagé s'il n'est pas utilisé correctement.

- Insérez l'embout pour le siège de l'aiguille complètement dans le tournevis dynamométrique.
- Insérez le tournevis dynamométrique avec l'embout pour le siège de l'aiguille droit dans le siège de l'aiguille.
- Ne poussez et ne tirez pas sur l'outil avec une force inutile.
- 5. Serrez tous les raccords sur le trajet d'écoulement à l'aide du tournevis dynamométrique muni de l'embout Viper.
- Exécutez le script D02 Tester le serrage du système (ce qui devrait être testé = l'échantillonneur) si la fuite continue.
- Des gouttelettes continuent à apparaître sur la vanne d'injection : Remplacez l'unité d'aiguille et le siège de l'aiguille. Reportez-vous à la section Remplacement du siège de l'aiguille et de l'unité de l'aiguille (▶ page 254).

7.15 Entretien du compartiment à colonnes

7.15.1 Remplacement de colonnes



ATTENTION—Colonnes chaudes

Les colonnes à l'intérieur de l'appareil peuvent devenir chaudes pendant l'utilisation. Ne touchez pas les surfaces chaudes pour éviter de vous brûler.

Laissez les colonnes refroidir avant de commencer les procédures de remplacement ou de maintenance.

À quel moment exécuter cette opération ?

Pour changer entre des colonnes

Éléments nécessaires

Colonne, selon les besoins de l'application

Préparatifs

1. Arrêtez le débit de la pompe.

Procédez comme suit

- 1. Retirez les capillaires de l'entrée et de la sortie de la colonne.
- 2. Retirez la colonne à remplacer de l'attache pour colonne. Pour retirer la bague de la partie inférieure de l'attache pour colonne, appuyez légèrement sur la bague dans la direction des flèches indiquées sur la bague et tirez-la vers vous.



Illustration 195: Retrait de la bague

- 3. Connectez les capillaires à l'entrée et à la sortie de la colonne de remplacement.
- 4. Placez la colonne dans la partie inférieure de l'attache pour colonne et réinstallez la bague.

5. Acheminez les capillaires par l'un des guides pour capillaires illustrés dans la figure.



Illustration 196: Guides pour capillaires

- Pour configurer les propriétés de la colonne, exécutez les scripts : A03 – Définir le type de colonne de séparation et A04 – Définir les caractéristiques de la colonne de séparation (et A05 – Définir les caractéristiques de la colonne de piégeage, le cas échéant).
- 7. Démarrez le débit de la pompe.

7.15.2 Inspection des vannes de commutation pour colonne à la recherche de fuites

La vanne Vanquish et la vanne VICI sont chacune munies d'un trou d'évacuation.



Illustration 197: Trous d'évacuation dans la vanne Vanquish (à gauche) et sur la vanne VICI (à droite)

À quel moment exécuter cette opération ?

- Des gouttelettes de liquide apparaissent au trou d'évacuation
- Du liquide s'est accumulé dans le bac de fuite

Procédez comme suit

- 1. Comme les fuites surviennent généralement aux raccords, inspectez visuellement tous les raccords du trajet d'écoulement.
- 2. Retirez puis réinstallez les capillaires.
- 3. Serrez tous les raccords sur le trajet d'écoulement.
- 4. Inspectez les vannes visuellement pour voir si elles fuient.
- Exécutez le script : D02 Tester le serrage du système (ce qui devrait être testé = l'échantillonneur). Ce script vérifie aussi si les vannes de commutation pour colonne présentent des fuites.
- S'il y a toujours une fuite, remplacez le joint de stator et/ou de rotor de la vanne Vanquish ou de la vanne VICI, en fonction de la vanne. Pour la vanne Vanquish, reportez-vous à la section Remplacement de la vanne de commutation Vanquish (▶ page 305).

Pour le joint du stator et/ou rotor de la vanne VICI, reportez-vous à la section Remplacement du joint du stator et/ou rotor de la vanne de commutation VICI () page 311).

7.15.3 Remplacement de la vanne de commutation Vanquish

À quel moment exécuter cette opération ?

- La vanne de commutation est défectueuse
- Pour changer entre différents types de vanne

Éléments nécessaires

- Vanne de commutation Vanquish, selon les exigences de l'application
- 4 vis

Outils nécessaires

Tournevis type Torx T10

Préparatifs

- 1. Arrêtez le débit de la pompe.
- 2. Déconnectez toutes les conduites de liquide raccordées à la vanne de commutation.
- 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 de remplacement.

Retrait de la vanne de commutation Vanquish

1. Desserrez les 4 vis de la vanne de commutation. Conservez les vis pour l'installation de la vanne de commutation de remplacement.



Illustration 198: 4 vis sur la vanne de commutation

 Retirez avec précaution la vanne de commutation du boîtier. Pour faciliter le retrait, vissez un bouchon de raccordement dans un orifice de vanne.





3. Déconnectez le câble du raccord situé à l'arrière de la vanne de commutation.



Illustration 200: Déconnexion du câble du raccord

Si aucune vanne n'est installée après le remplacement

- 1. Scellez le connecteur de câbles à l'aide de ruban adhésif étanche.



- 2. Enfoncez le câble dans le boîtier.
- 3. Placez le couvercle sur l'ouverture de la vanne.
- 4. Serrez les 4 vis du couvercle.

Installation de la vanne de commutation Vanquish

1. Connectez la vanne de commutation de remplacement au câble.



Illustration 202: Connexion du câble

2. Enfoncez le câble dans le boîtier.



3. Enfoncez délicatement la vanne de commutation dans le boîtier.

Illustration 203: Installation de la vanne de commutation Vanquish

- 4. Vérifiez que la vanne de commutation est installée dans la bonne direction : l'orifice n° 3 doit être situé en haut.
- 5. À l'aide d'un tournevis, serrez les 4 vis de la vanne de commutation.
- 6. Retirez les bouchons de raccordement qui couvrent les orifices des vannes.
- 7. Connectez les conduites de liquide, selon les besoins de l'application.
- 8. Mettez le compartiment à colonnes sous tension.
- 9. Inspectez les vannes visuellement pour voir si elles fuient.
- Exécutez le script : D02 Tester le serrage du système (ce qui devrait être testé = l'échantillonneur). Ce script vérifie aussi si les vannes de commutation pour colonne présentent des fuites.

7.15.4 Remplacement de la vanne de commutation VICI

Vous ne pouvez remplacer la vanne VICI qu'aux emplacements où une commande de vanne VICI est installée. Pour installer la commande de vanne VICI, prenez contact avec l'assistance technique de Thermo Fisher Scientific locale. Pour une vue d'ensemble de la vanne de commutation VICI, reportez-vous à la section Vue d'ensemble de la vanne de commutation pour colonne (▶ page 54).

Prérequis

La commande de la vanne VICI et le cadre d'adaptateur sont installés (effectué par l'ingénieur d'entretien).

À quel moment exécuter cette opération ?

- La vanne de commutation est défectueuse
- Pour changer entre différents types de vanne

Éléments nécessaires

• Vanne de commutation VICI selon les exigences de l'application

Outils nécessaires

Tournevis type Torx T10

Préparatifs

- 1. Arrêtez le débit de la pompe.
- 2. Déconnectez toutes les conduites de liquide raccordées à la vanne de commutation VICI.
- Ne retirez pas encore les bouchons de raccordement qui couvrent les orifices des vannes sur la vanne de commutation de remplacement.
- Pour retirer la capsule de vanne installée de la bride de la vanne, vissez un bouchon de raccordement dans un orifice de vanne et tirer fermement sur la capsule de vanne à l'aide du bouchon de raccordement.

Si aucune vanne n'est installée après le remplacement

- 1. Desserrez les 4 vis du cadre de l'adaptateur.
- 2. Retirez le cadre de l'adaptateur.
- 3. Placez le couvercle sur l'ouverture de la vanne.
- 4. Serrez les 4 vis du couvercle.

Installation de la vanne de commutation VICI

 Alignez les deux broches situées à l'arrière de la bague de verrouillage de la vanne avec les ouvertures correspondantes sur la bride de la vanne.



Illustration 204: Alignement de la bague de verrouillage et de la bride

- 2. Pressez la capsule de la vanne sur la bride de la vanne.
- Pour aligner l'axe avec la capsule, appuyez sur le bouton VALVE LEFT ou le bouton VALVE RIGHT sur le panneau de commande et maintenez-le enfoncé, en fonction de la position de la vanne. La fonction de vanne de captage s'active et vous pouvez relâcher le bouton.
- 4. Quand l'axe est aligné, la capsule de vanne s'enclenche dans la bride de la vanne.
- 5. Retirez les bouchons de raccordement qui couvrent les orifices des vannes.
- 6. Connectez les conduites de liquide, selon les besoins de l'application.
- 7. Configurez la vanne de commutation VICI dans le système de gestion de données chromatographiques.
- 8. Inspectez les vannes visuellement pour voir si elles fuient.
- Exécutez le script : D02 Tester le serrage du système (ce qui devrait être testé = l'échantillonneur). Ce script vérifie aussi si les vannes de commutation pour colonne présentent des fuites.

7.15.5 Remplacement du joint du stator et/ou rotor de la vanne de commutation VICI

À quel moment exécuter cette opération ?

- Fuite de la vanne de commutation causée par un joint de stator et/ ou rotor endommagé
- Encrassement du joint du rotor

Éléments nécessaires

- Joint de remplacement du stator et/ou rotor de la vanne
- Connecteur Viper ou bouchons de raccordement en plastique

Outils nécessaires

Clé hexagonale, taille 9/64 po

Préparatifs

- 1. Arrêtez le débit de la pompe.
- 2. Déconnectez toutes les conduites de liquide raccordées à la vanne de commutation VICI.
- 3. Pour retirer la capsule de vanne installée de la bride de la vanne, vissez un connecteur Viper dans la capsule de vanne ou utilisez des bouchons de raccordement en plastique.
- 4. Tirez fermement sur la capsule de vanne à l'aide du connecteur Viper ou des bouchons de raccordement en plastique.

Procédez comme suit

1. Retirez les 2 vis du stator à l'aide de la clé hexagonale. Autrement, desserrez les vis l'une après l'autre jusqu'à ce que vous pouvez les retirer.



Illustration 205: Vis du stator

2. Retirez le stator du corps de la vanne. Pour éviter d'endommager la surface de scellage, tournez le stator pour qu'il repose sur la surface externe.



Illustration 206: vanne de commutation VICI

N°	Description
1	Stator
2	Corps de la vanne

3. Retirez avec précaution le joint du rotor manuellement.

AVIS

Pour éviter de rayer les surfaces de scellage, n'utilisez pas des outils pointus ou coupants pour retirer ou installer le joint. Même des petites rayures peuvent avoir des répercussions négatives sur les performances de la vanne.

- Vérifiez s'il y a des rayures sur les surfaces de scellage du joint du rotor et du stator. Si des rayures sont visibles, remplacez le joint du rotor et/ou le stator.
- 5. Insérez le joint de rotor de remplacement :
 - Respectez l'orientation du joint. La surface de scellage où sont gravés les passages de flux doit être à l'extérieur. Le profil est asymétrique pour empêcher l'installation incorrecte du joint du rotor.
 - Évitez soigneusement toute contamination des éléments de la vanne. Même de petites particules peuvent l'endommager et résulter en une mauvaise performance de scellage.
- 6. Réinstallez le stator et serrez les deux vis en alternance, environ un quart de tour chaque fois, jusqu'à ce qu'elles soient serrées.

NOTE Évitez de les serrer de façon excessive. Les vis maintiennent l'assemblage ensemble et ne touchent pas la force de scellage qui est réglée automatiquement quand les vis ferment le stator contre le corps de la vanne.

- Réinstallez la capsule de vanne tel que décrit dans la section respective dans Remplacement de la vanne de commutation VICI (> page 308).
- 8. Inspectez les vannes visuellement pour voir si elles fuient.
- Exécutez le script : D02 Tester le serrage du système (ce qui devrait être testé = l'échantillonneur). Ce script vérifie aussi si les vannes de commutation pour colonne présentent des fuites.
- Si vous constatez toujours une fuite, remplacez la vanne. Reportezvous à la section Remplacement de la vanne de commutation VICI (> page 308).

7.15.6 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.
- 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 92).

7.16 Contrôleur du système et maintenance du dispositif d'affichage

La présente section décrit comment entretenir le contrôleur du système et le dispositif d'affichage.

7.16.1 Remplacement du contrôleur du système

- 1. Retirez le contrôleur du système. Reportez-vous à la section Retrait du contrôleur du système (▶ page 314).
- Installez le contrôleur du système. Reportez-vous à la section Installation du contrôleur du système (> page 315).
- 3. Connectez le contrôleur du système. Reportez-vous à la section Connexion du contrôleur du système (▶ page 318).
- Si vous avez remplacé le contrôleur du système : reconfigurez votre système dans l'interface utilisateur Vanquish. Pour obtenir plus d'informations sur les scripts de configuration, reportez-vous à la section Préparation du système pour l'utilisation initiale (> page 139).

7.16.1.1 Retrait du contrôleur du système

Procédez comme suit

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

- Pour déplacer légèrement le socle de système et mieux accéder aux pièces du système, déverrouillez le socle du système à l'aide des outils de déverrouillage du socle. Reportez-vous à la section Déverrouillage du socle du système (> page 73).
- 2. Placez le système empilé dans une position de travail appropriée.
- 3. Débranchez les câbles de signal et d'alimentation du contrôleur du système.

4. Appuyez simultanément sur les deux agrafes-ressorts et retirez fermement le contrôleur du système.



Illustration 207: Pression sur les agrafes-ressorts du côté gauche et du côté droit du contrôleur du système

7.16.1.2 Installation du contrôleur du système

Éléments nécessaires

Bouton d'alimentation VSC

Outils nécessaires

- Paire de pinces
- Facultatif : Tournevis type Torx T10

Procédez comme suit

La fente prévue pour le contrôleur du système se trouve à l'arrière du socle du système.

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

- Pour déplacer légèrement le socle de système et mieux accéder aux pièces du système, déverrouillez le socle du système à l'aide des outils de déverrouillage du socle. Reportez-vous à la section Déverrouillage du socle du système (▶ page 73).
- 2. Placez le système empilé dans une position de travail appropriée.
- 3. Insérez la carte SD en respectant les consignes suivantes :
 - Orientation correcte de l'encoche (reportez-vous à l'image)
 - Insérer la carte SD droit dans le système de fixation intérieur avant de l'enfoncer complètement dans la fente.



Illustration 208: Insertion de la carte SD

 À l'arrière du système empilé, soulevez légèrement le contrôleur du système et enfoncez-le fermement dans le socle du système jusqu'à ce qu'il s'enclenche en position.



Illustration 209: Insertion du contrôleur du système

5. Si le contrôleur du système ne s'enclenche pas en position même si vous l'enfoncez fermement : Retirez le support de montage en desserrant les 3 vis avec le tournevis.



Illustration 210: Retrait du support de montage

N°	Description
1	Support de montage
2	Vis

AVIS—Dommages au contrôleur du système

Sans support de montage, le contrôleur du système peut tomber et être endommagé quand vous soulevez le système. Cela pourrait endommager le contrôleur du système.

- Connectez les câbles d'interface. Reportez-vous à la section Connexion du contrôleur du système (▶ page 318).
- Si vous avez remplacé le contrôleur du système : reconfigurez votre système dans l'interface utilisateur Vanquish. Pour obtenir plus d'informations sur les scripts de configuration, reportez-vous à la section Préparation du système pour l'utilisation initiale (> page 139).

7.16.1.3 Connexion du contrôleur du système

Les connecteurs et emplacements suivants sont fournis sur le panneau court du contrôleur du système :



Illustration 211: Connecteur sur le panneau court

N°	Description
1	Plaque signalétique
	Indique le numero de serie et la tension nominale.
2	Connecteur Vanquish System Controller (VSC)
	Permet la connexion au socle du système.
3	Port System Interlink
	Permet le raccordement au socle du système et à un module dans le système HPLC
4	Port Universal Serial Bus (USB) (connecteur de type "A")
	Désactivé – Ne pas utiliser
5	Port Local Area Network (LAN)
	Permet la connexion au réseau du site.
6	Interrupteur d'alimentation principal (marche/arrêt)
7	Port d'entrée d'alimentation

Les connecteurs et emplacements suivants sont fournis sur le panneau long du contrôleur du système :



Illustration 212: Connecteurs sur le panneau long

N°	Description
1	Port de sortie d'alimentation du dispositif d'affichage
2	Port de données du dispositif d'affichage
3	Ports System Interlink Désactivés
4	Port USB (connecteur de type "A") Désactivé – Ne pas utiliser
5	Étiquette indiquant le numéro de référence, le nom du module et le numéro de révision.
6	Emplacement pour carte mémoire flash (SD)

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 le système.

Éléments nécessaires

- Cordon d'alimentation
- Câble VSC court (longueur : 8 cm)
- Câble System Interlink (longueur : 50 cm)
- Si vous souhaitez accéder au contrôleur du système à partir du réseau de votre site : câble LAN (longueur : 5 m)

Procédez comme suit

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.
- À l'aide du câble VSC court, connectez le port VSC du socle du système au port VSC du contrôleur du système.



Illustration 213: Connexion des câbles VSC et System Interlink

AVIS—Endommagement du module Vanquish

Les ports **System Interlink** et **LAN** utilisent le même type de connecteurs. Le raccordement d'un câble LAN, par exemple, au port **System Interlink** d'un module Vanquish, peut endommager ce module et le contrôleur du système.

Branchez uniquement les câbles sur les ports correspondants.

- 5. À l'aide d'un câble System Interlink, connectez le port System Interlink du socle du système à un port System Interlink sur le panneau court du contrôleur du système.
- 6. Connectez l'autre port System Interlink du panneau court du contrôleur du système au port System Interlink de la pompe.
- 7. Branchez un câble LAN sur le port LAN du contrôleur du système.
- 8. Branchez l'autre connecteur du câble LAN sur le port LAN de carte réseau de l'ordinateur sur lequel est installé le système de gestion de données chromatographiques.
- 9. Replacez le système empilé dans la position d'utilisation.
- Verrouillez le socle du système à l'aide des outils de verrouillage (reportez-vous à la section Verrouillage du socle du système (> page 119)).

7.16.2 Remplacement de la carte SD

NOTE Pour être sûr que la carte SD est compatible avec le contrôleur du système, utilisez uniquement les cartes SD commandées auprès de Thermo Fisher Scientific.

À quel moment exécuter cette opération ?

Un message s'affiche sur la Vanquish User Interface pour indiquer que la carte SD doit être remplacée.

Matériel nécessaire

- Nouvelle carte SD
- En option : crayon ou tournevis plat

Préparatifs

- 1. Redémarrez le contrôleur du système et attendez que la LED d'état du contrôleur s'allume.
- 2. Vérifiez si le message réapparaît. Si c'est le cas, passez à l'étape suivante.
- 3. Arrêtez le contrôleur du système au moyen du bouton marche/arrêt du contrôleur du système.

Procédez comme suit

- Pour déplacer légèrement le socle de système et mieux accéder aux pièces du système, déverrouillez le socle du système à l'aide des outils de déverrouillage du socle. Reportez-vous à la section Déverrouillage du socle du système (> page 73).
- 2. Placez le système empilé dans une position de travail appropriée.
- 3. Appuyez légèrement sur la carte SD jusqu'à ce qu'elle se libère. Si nécessaire, utilisez un crayon ou un tournevis plat.



Illustration 214: Retrait de la carte SD

- 4. Retirez l'ancienne carte SD.
- 5. Insérez la carte SD en respectant les consignes suivantes :
 - Orientation correcte de l'encoche (reportez-vous à l'image)
 - Insérer la carte SD droit dans le système de fixation intérieur avant de l'enfoncer complètement dans la fente.





6. Replacez le système empilé dans la position d'utilisation.

 Verrouillez le socle du système à l'aide des outils de verrouillage (reportez-vous à la section Verrouillage du socle du système (> page 119)).

7.16.3 Remplacement du dispositif d'affichage

- 1. Désinstallez l'ancien dispositif d'affichage. Reportez-vous à la section Désinstallation du dispositif d'affichage (▶ page 323).
- 2. Installez le nouveau dispositif d'affichage. Reportez-vous à la section Installation du dispositif d'affichage (▶ page 99).
- 3. Connectez les câbles du dispositif d'affichage. Reportez-vous à la section Connexion des câbles du dispositif d'affichage (▶ page 111).

7.16.3.1 Désinstallation du dispositif d'affichage

Outils nécessaires

Tournevis type Torx T10

Procédez comme suit

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

- Pour déplacer légèrement le socle de système et mieux accéder aux pièces du système, déverrouillez le socle du système à l'aide des outils de déverrouillage du socle. Reportez-vous à la section Déverrouillage du socle du système (> page 73).
- 2. Placez le système empilé dans une position de travail appropriée.
- 3. Débranchez les deux connecteurs du dispositif d'affichage du contrôleur du système.
- 4. Retirez le cache du câble.
- 5. Retirez le câble du dispositif d'affichage du rail et des serre-câbles du boîtier.

6. Tout en maintenant le bloc de fixation avec la main droite, desserrez la vis d'un des trous marqués d'un cercle de l'extérieur avec la main gauche.



Illustration 216: Desserrage de la vis maintenant en place le bloc de fixation

- 7. Retirez le bloc de fixation.
- 8. Desserrez les 2 vis dans les deux trous supérieurs.



Illustration 217: Desserrage des vis

9. Poussez l'assemblage du dispositif d'affichage vers l'arrière du système et retirez-le du système empilé.


10. Fixez les 2 vis supérieures et fixez le bloc de fixation avec une vis.

Illustration 218: Vis à fixer sur l'assemblage du dispositif d'affichage

7.17 Remplacement du bac à solvants

Éléments nécessaires

- Bac à solvant de remplacement
- Serviette en papier non pelucheuse

Outils nécessaires

Tournevis type Torx T10

Préparatifs

- 1. Arrêtez le système au moyen du bouton d'alimentation du système (bouton du bas sur le socle du système).
- Arrêtez le contrôleur du système au moyen de son bouton d'alimentation électrique (bouton du haut sur le socle du système).
- 3. Videz le système de rinçage des joints :
 - a) Dévissez le bouchon du réservoir de liquide de rinçage des joints.
 - b) Retirez la conduite de rinçage des joints en même temps que le bouchon du réservoir de liquide de rinçage des joints.
 - c) Pour vider la conduite de rinçage, soulevez le levier de la pompe de rinçage des joints. Relâchez le levier lorsque la conduite est vide.
- 4. Déconnectez et videz les conduites de solvant :
 - a) Dévissez les bouchons des réservoirs de solvant.
 - b) Retirez en même temps les bouchons et les conduites de solvant des réservoirs.
 - c) Protégez les filtres de conduite de solvant de toute contamination de manière appropriée.
 - d) Déconnectez les conduites de solvant des orifices d'admission la pompe et placez l'extrémité libre des conduites de solvant dans le bac de fuite pour vider les conduites.
 - e) Fermez les orifices d'aspiration de solvant de la pompe avec des capuchons appropriés.
- 5. Absorbez tout liquide des conduites de liquides avec une serviette en papier non pelucheuse.
- 6. Retirez les réservoirs de liquides du bac à solvants.

Procédez comme suit

- 1. Desserrez les vis qui fixent le support de montage au module du système sous le bac à solvants, à l'aide du tournevis.
- 2. Retirez le support en tôle.

NOTE Gardez le support et les vis. Vous pourriez en avoir besoin pour un transport ultérieur.



Illustration 219: Retrait du support en tôle (ici : passeur d'échantillon)

- 3. Soulevez la partie avant du bac à solvants et retirez-le en tirant vers l'avant.
- 4. Tirez les conduites de liquide en dehors de la gaine de tubulure du bac à solvants.
- Retirez l'insertion de bouteille du bac à solvants (reportez-vous à la section Retrait de l'insertion de bouteille (▶ page 329)) si disponible.
- 6. Faites passer les conduites de liquide à travers la gaine de tubulure du nouveau bac à solvants.
- Placez le nouveau bac à solvants sur le module du système du haut, environ 5 cm avant la fin des rails. Poussez le bac à solvants vers l'arrière jusqu'à ce qu'il s'enclenche en position.
- Positionnez l'insertion de bouteille dans le bac à solvants (reportezvous à la section Installation de l'insertion de bouteille (> page 330)) si disponible.
- 9. Placez les réservoirs à liquide sur le bac à solvants.
- 10. Connectez les conduites de liquide aux réservoirs de liquides.
- 11. Raccordez les conduites de solvant aux orifices d'admission de la pompe.
- 12. Mettez le contrôleur du système sous tension au moyen de son bouton d'alimentation électrique (bouton du haut sur le socle du système).

- 13. Mettez le système sous tension au moyen du bouton d'alimentation du système (bouton du bas sur le socle du système).
- 14. Préparez le système pour l'utilisation initiale. Reportez-vous à la section Préparation du système pour l'utilisation initiale (▶ page 139).

7.18 Remplacement de l'insertion de bouteille

7.18.1 Retrait de l'insertion de bouteille

Éléments nécessaires

Serviette en papier non pelucheuse

Procédez comme suit

- 1. Arrêtez le système au moyen du bouton d'alimentation électrique (bouton du bas sur le socle du système).
- Arrêtez le contrôleur du système au moyen de son bouton d'alimentation électrique (bouton du haut sur le socle du système).
- 3. Déconnectez les conduites de rinçage des joints :
 - a) Dévissez le bouchon du réservoir de liquide de rinçage des joints.
 - b) Retirez la conduite de rinçage des joints en même temps que le bouchon du réservoir de liquide de rinçage des joints.
- 4. Déconnectez les conduites de solvant :
 - a) Dévissez les bouchons des réservoirs de solvant.
 - b) Retirez en même temps les bouchons et les conduites de solvant des réservoirs.
 - c) Protégez les filtres de conduite de solvant de toute contamination de manière appropriée.
- 5. Absorbez tout liquide des conduites de liquides avec une serviette en papier non pelucheuse.
- 6. Retirez les réservoirs de liquides du bac à solvants.
- 7. Poussez fermement les ailes de l'insertion de bouteille au-dessus du rail du bac à solvants et retirez l'insertion de bouteille.



Illustration 220: Retrait de l'insertion de bouteille

N°	Description
1	Aile

7.18.2 Installation de l'insertion de bouteille

Éléments nécessaires

Insertion de bouteille

Procédez comme suit

- 1. Insérez l'aile arrière et gauche dans l'espace en dessous des rails.
- 2. Poussez fermement l'aile droite en dessous des rails.



Illustration 221: Insertion de l'insertion de bouteille

- 3. Placez les réservoirs à liquide dans l'insertion de bouteille du bac à solvants.
- 4. Connectez les conduites de liquide aux réservoirs de liquides.

- Mettez le système et le contrôleur du système sous tension au moyen du bouton d'alimentation situé à gauche du socle du système.
- Préparez le système pour l'utilisation initiale. Reportez-vous à la section Préparation du système pour l'utilisation initiale (> page 139).

7.19 Remplacement des fusibles de l'alimentation principale et des portes

7.19.1 Remplacement des fusibles de l'alimentation principale

À quel moment exécuter cette opération ?

Fusibles grillés

Éléments nécessaires

Fusibles (2 fusibles, 5 AT, 230 V AC, fusion lente, 5 \times 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 222: 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.

7.19.2 Remplacement des portes

À quel moment exécuter cette opération ?

Porte endommagée

NOTE Les procédures de maintenance ne nécessitent pas de retirer les portes. Si cela est nécessaire pour une raison ou une procédure en particulier, suivez les étapes de cette section.

Éléments nécessaires

Porte de remplacement

Préparatifs

Si la porte à remplacer est située directement en dessous du bac à solvants, procédez comme suit :

- 1. Retirez tous les réservoirs du bac à solvants.
- Retirez le support en tôle qui fixe le bac à solvants au passeur d'échantillon, le cas échéant. Pour les instructions, reportez-vous à la section Installation du compartiment à colonnes dans le système empilé (> page 86).

Si le dispositif d'affichage est installé près de la porte à remplacer, envisagez de desserrer le bloc de fixation du dispositif d'affichage pour pousser le dispositif vers l'arrière du système.

Pour obtenir des informations sur comment desserrer le bloc de fixation du dispositif d'affichage, reportez-vous à la section Désinstallation du dispositif d'affichage (▶ page 323).

Procédez comme suit

AVIS

Pour éviter d'endommager les charnières de la porte, soyez prudent lorsque vous exécutez la procédure suivante et ne forcez pas.

- 1. Si la porte est placée directement en dessous du bac à solvants, soulevez légèrement le rebord avant du bac.
- 2. Pour retirer une porte, poussez-la vers le haut pendant son ouverture.

Placez la porte de façon à aligner les deux charnières du boîtier sur les rainures de la porte. Vous pouvez décrocher la porte seulement si les charnières sont dans les rainures.



Illustration 223: Décrochage d'une porte

N°	Description
1	Charnières du boîtier
2	Rainure de réception sur la porte

- 3. Inclinez légèrement la porte vers l'extérieur du boîtier, puis retirezla.
- 4. Pour installer la porte, alignez-la sur les charnières du boîtier. Ne coincez pas les tubes ou les capillaires entre la porte et le boîtier.
- 5. Insérez les charnières dans la rainure en exerçant une pression vers le haut et en tournant légèrement la porte.
- Poussez la porte vers le bas pour la verrouiller en place. Vous pouvez fermer la porte seulement si elle est correctement installée.

7.20 Remplacement du module à glissières

Vous pouvez retirer le module à glissières du boîtier d'un module en vue du transport ou de l'expédition. Le boîtier reste dans le système empilé. Pour retourner un module défectueux à l'usine, installez le module à glissières dans le boîtier du module de remplacement.

7.20.1 Préparation du retrait de la pompe

Préparatifs généraux

- 1. Exécutez le script : **B05 Arrêt (de longue durée)** avec le paramètre suivant : **Module à retirer = pompe**.
- Pour déplacer légèrement le socle de système et mieux accéder aux pièces du système, déverrouillez le socle du système à l'aide des outils de déverrouillage du socle. Reportez-vous à la section Déverrouillage du socle du système (▶ page 73).
- 3. Placez le système empilé dans une position de travail appropriée.

Arrêt

- Mettez *tous* les modules du système hors tension au moyen de leurs interrupteurs principaux d'alimentation. Il ne suffit pas d'appuyer sur le bouton d'alimentation du système pour couper complètement l'alimentation électrique des appareils.
- 2. Débranchez le cordon d'alimentation.
- 3. Débranchez tous les câbles de signal vers d'autres appareils.

Raccordements fluidiques

- 1. Retirez tous les raccordements fluidiques vers d'autres modules du système :
 - Conduite d'admission de la pompe de rinçage des joints du tube péristaltique de la pompe de rinçage des joints
 - Conduites de rinçage de l'aiguille (étiquetées W et S) des orifices du dégazeur
 - Capillaire vers le passeur d'échantillon de l'orifice de refoulement de la pompe
 - Conduites de solvant des orifices d'admission de la pompe

Les raccordements fluidiques et raccords de capillaires sont extrêmement sensibles à la contamination. La poussière et les débris peuvent contaminer ces raccordements.

Installez toujours des capuchons sur les capillaires et des bouchons sur les raccordements fluidiques ouverts pour éviter toute contamination.

- 2. Fermez les raccordements fluidiques ouverts au moyen de bouchons appropriés, par exemple les bouchons du kit d'expédition.
- 3. Vérifiez le bac de fuite et l'orifice d'évacuation. S'il y a du liquide, absorbez-le avec une serviette en papier.

7.20.2 Retrait du module à glissières de la pompe



ATTENTION—Charge lourde, instrument encombrant

Les modules du système sont trop lourds ou encombrants pour être soulevés en toute sécurité par une seule personne. Afin d'éviter toute blessure corporelle ou détérioration des modules du système, veuillez respecter les consignes suivantes :

- Le maniement physique des modules, notamment son soulèvement ou son déplacement, requiert la présence d'au moins deux personnes.
- La présence de plusieurs personnes est requise en particulier pour placer les modules dans le système empilé ou pour les en retirer.

Outils nécessaires

Tournevis type Torx T20

Préparatifs

Reportez-vous à la section Préparation du retrait de la pompe (> page 336).

Procédez comme suit

1. Desserrez les quatre vis imperdables situées sur le devant gauche et droit de l'appareil.



Illustration 224: Vis imperdables du module à glissières de la pompe

- Poussez tous les tubes et capillaires présents dans la gaine de tubulure des modules du système Vanquish, dans la gaine de tubulure.
- 3. Saisissez le module à glissières par le bac de fuite, puis retirez le module du boîtier d'environ 10 cm.

Le module à glissières peut tomber si vous le retirez trop du boîtier.

Tirez sur le module à glissières juste assez pour pouvoir le saisir par les deux côtés et par le bas.

- 4. Retirez le module à glissières du boîtier. La procédure suivante requiert la présence d'au moins deux personnes :
 - a) Saisissez le module à glissières par le bas des deux côtés.
 - b) Retirez le module à glissières des rails en le tirant vers l'avant.
 - c) Placez le module à glissières sur une surface propre et stable.

7.20.3 Préparation du retrait du passeur d'échantillon

Préparatifs généraux

- 1. Retirez les portoirs d'échantillons et les récipients d'échantillonnage du compartiment à échantillons.
- Exécutez le script : B05 Arrêt (de longue durée) avec le paramètre suivant : Module à retirer = échantillonneur.
- 3. Si ce n'est pas déjà fait, verrouillez l'unité d'aiguille (reportez-vous à la section Verrouillage de l'unité d'aiguille (▶ page 251)).

- Pour déplacer légèrement le socle de système et mieux accéder aux pièces du système, déverrouillez le socle du système à l'aide des outils de déverrouillage du socle. Reportez-vous à la section Déverrouillage du socle du système (▶ page 73).
- 5. Placez le système empilé dans une position de travail appropriée.

Arrêt

- Mettez tous les modules du système hors tension au moyen de leurs interrupteurs principaux d'alimentation. Il ne suffit pas d'appuyer sur le bouton d'alimentation du système pour couper complètement l'alimentation électrique des appareils.
- 2. Débranchez le cordon d'alimentation.
- 3. Débranchez tous les câbles de signal vers d'autres appareils.

Raccordements fluidiques

- 1. Retirez tous les raccordements fluidiques vers d'autres modules du système et aux réservoirs :
 - Conduite du réservoir de liquide de rinçage des joints et conduite d'admission de la pompe de rinçage des joints. Raccordez les conduites qui restent connectées à la tête du dispositif de mesure avec un raccord de tube.
 - Conduites de rinçage de l'aiguille
 - Capillaire depuis la pompe
 - Capillaire vers le compartiment à colonnes, si installé

AVIS

Les raccordements fluidiques et raccords de capillaires sont extrêmement sensibles à la contamination. La poussière et les débris peuvent contaminer ces raccordements.

Installez toujours des capuchons sur les capillaires et des bouchons sur les raccordements fluidiques ouverts pour éviter toute contamination.

Nettoyage

- En présence de condensation ou si des échantillons ont été renversés dans le compartiment à échantillons, nettoyez et décontaminez le compartiment à échantillons avant de poursuivre.
- 2. Vérifiez si les bacs de fuite contiennent des liquides provenant de fuites. Si c'est le cas, absorbez-les avec une serviette en papier.

7.20.4 Retrait du module à glissières du passeur d'échantillon



ATTENTION—Charge lourde, instrument encombrant

Les modules du système sont trop lourds ou encombrants pour être soulevés en toute sécurité par une seule personne. Afin d'éviter toute blessure corporelle ou détérioration des modules du système, veuillez respecter les consignes suivantes :

- Le maniement physique des modules, notamment son soulèvement ou son déplacement, requiert la présence d'au moins deux personnes.
- La présence de plusieurs personnes est requise en particulier pour placer les modules dans le système empilé ou pour les en retirer.

Outils nécessaires

Tournevis type Torx T20

Préparatifs

Reportez-vous à la section Préparation du retrait du passeur d'échantillon () page 338).

Retrait du module à glissières

1. Desserrez les quatre vis imperdables situées sur le devant gauche et droite de l'appareil.



Illustration 225: Vis imperdables du module à glissières du passeur d'échantillon

- 2. Poussez tous les tubes et capillaires présents dans la gaine de tubulure des modules du système Vanquish, dans la gaine de tubulure.
- 3. Saisissez le module à glissières par le cadre central où se trouve le verrou de la porte, puis retirez le module du boîtier d'environ 10 cm.

Le module à glissières peut tomber si vous le retirez trop du boîtier.

Tirez sur le module à glissières juste assez pour pouvoir le saisir par les deux côtés et par le bas.

- 4. Retirez le module à glissières du boîtier. La procédure suivante requiert la présence d'au moins deux personnes :
 - a) Saisissez le module à glissières par le bas des deux côtés.
 - b) Retirez le module à glissières des rails en le tirant vers l'avant.
 - c) Placez le module à glissières sur une surface propre et stable.

7.20.5 Préparation du retrait du compartiment à colonnes

Préparatifs généraux

- 1. Exécutez le script : **B05 Arrêt (de longue durée)** avec le paramètre suivant : **Module à supprimer = Compartiment à colonnes**.
- Pour déplacer légèrement le socle de système et mieux accéder aux pièces du système, déverrouillez le socle du système à l'aide des outils de déverrouillage du socle. Reportez-vous à la section Déverrouillage du socle du système (▶ page 73).
- 3. Placez le système empilé dans une position de travail appropriée.

Arrêt

- 1. Mettez *tous* les modules du système hors tension au moyen de leurs interrupteurs principaux d'alimentation. Il ne suffit pas d'appuyer sur le bouton d'alimentation du système pour couper complètement l'alimentation électrique des appareils.
- 2. Débranchez le cordon d'alimentation.
- 3. Débranchez tous les câbles de signal vers d'autres appareils.

Raccordements fluidiques

- 1. Retirez tous les raccordements fluidiques vers d'autres modules du système.
- 2. Couvrez les orifices ouverts des vannes de commutation avec de bouchons de raccordement.

Les raccordements fluidiques et raccords de capillaires sont extrêmement sensibles à la contamination. La poussière et les débris peuvent contaminer ces raccordements.

Installez toujours des capuchons sur les capillaires et des bouchons sur les raccordements fluidiques ouverts pour éviter toute contamination.

Accessoires en option

- Retirez la colonne en suivant les instructions de la section Installation des colonnes dans le compartiment à colonnes (> page 94) dans le sens inverse.
- Retirez l'appareil de préchauffage, s'il est installé (voir Remplacement de l'appareil de préchauffage passif (> page 313)).
- Retirez les vannes de commutation (reportez-vous à la section Remplacement de la vanne de commutation Vanquish (▶ page 305) ou Remplacement de la vanne de commutation VICI (▶ page 308)).

Nettoyage

Vérifiez le bac de fuite et l'orifice d'évacuation. Si c'est le cas, absorbezles avec une serviette en papier.

7.20.6 Retrait du module à glissières du compartiment à colonnes

Outils nécessaires

Tournevis type Torx T20

Préparatifs

Reportez-vous à la section Préparation du retrait du compartiment à colonnes () page 341).

Retrait du module à glissières

1. Desserrez les quatre vis imperdables situées sur le devant gauche et droite de l'appareil.



Illustration 226: Vis imperdables du module à glissières du compartiment à colonnes

- 2. Poussez tous les tubes et capillaires présents dans la gaine de tubulure des modules du système Vanquish, dans la gaine de tubulure.
- 3. Saisissez le module à glissières par le bac de fuite, puis retirez le module du boîtier d'environ 10 cm.

Le module à glissières peut tomber si vous le retirez trop du boîtier.

Tirez sur le module à glissières juste assez pour pouvoir le saisir par les deux côtés et par le bas.

- 4. Retirez le module à glissières du boîtier. La procédure suivante requiert la présence d'au moins deux personnes :
 - a) Saisissez le module à glissières par le bas des deux côtés.
 - b) Retirez le module à glissières des rails en le tirant vers l'avant.
 - c) Placez le module à glissières sur une surface propre et stable.

7.20.7 Retour du module à glissières



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.

Préparatifs

Retirez le module à glissières que vous souhaitez retourner du boîtier. Voir l'information pertinente :

Retrait du module à glissières de la pompe () page 337)

Retrait du module à glissières du passeur d'échantillon () page 340)

Retrait du module à glissières du compartiment à colonnes () page 342)

Procédez comme suit

- Installez le module à glissières dans le boîtier de l'appareil de remplacement. Suivez les instructions relatives à l'insertion du module à glissières dans le boîtier et au serrage des vis imperdables de la section Installation du module à glissières (> page 345).
- 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.

7.20.8 Installation du module à glissières



ATTENTION—Charge lourde, instrument encombrant

Les modules du système sont trop lourds ou encombrants pour être soulevés en toute sécurité par une seule personne. Afin d'éviter toute blessure corporelle ou détérioration des modules du système, veuillez respecter les consignes suivantes :

- Le maniement physique des modules, notamment son soulèvement ou son déplacement, requiert la présence d'au moins deux personnes.
- La présence de plusieurs personnes est requise en particulier pour placer les modules dans le système empilé ou pour les en retirer.

Outils nécessaires

Tournevis type Torx T20

Préparatifs

1. Lorsque vous installez le module à glissières dans un boîtier du système empilé, vérifiez que ce boîtier est correctement placé dans la pile.

Procédez comme suit

- Poussez tous les tubes et capillaires présents dans la gaine de tubulure des modules du système Vanquish Neo, dans la gaine de tubulure. Dans le cas contraire, vous ne pourrez pas, à l'étape suivante, insérer correctement le module à glissières dans le boîtier.
- 2. Dans le module à glissières, vérifiez que vous pouvez déplacer les vis imperdables d'avant en arrière avec vos doigts. Si cela n'est pas le cas, vissez-les de manière à pouvoir le faire.
- 3. Insérez le module à glissières dans le boîtier. Les étapes suivantes nécessitent un effort d'équipe :
 - a) Saisissez le module à glissières par le bas des deux côtés.
 - b) Soulevez le module à glissières à hauteur du boîtier.
 - c) Placez le module à glissières dans le boîtier de telle sorte qu'il y soit inséré d'environ 25 cm.
 - d) Enfoncez le module à glissières sur les rails et dans le boîtier jusqu'à ce qu'il soit complètement inséré dans le boîtier.



Illustration 227: Insertion du module à glissières

- 4. Pour faciliter le vissage du module à glissières au boîtier, tirez et sortez le module d'environ 5 mm.
- Avec le tournevis, déplacez chaque vis imperdable d'avant en arrière tout en poussant vers l'intérieur jusqu'à ce que la vis glisse dans le filetage.
- 6. Réinsérez le module à glissières dans le boîtier aussi loin que possible.
- 7. Serrez manuellement les quatre vis imperdables du module à glissières de manière progressive et uniforme.

- Assurez-vous que les vis sont serrées. Tirez le module à glissières vers l'avant par le bac de fuite pour vérifier si les vis bougent. Si elles ne bougent pas, alors le module à glissières est correctement installé.
- Si les vis bougent, serrez-les davantage. Avec une clé dynamométrique, le couple recommandé est de 1,2 Nm.
- Pour configurer le module, suivez les étapes de la section Installation (> page 75).
- 9. Faites programmer le numéro de série du système dans le module du système par un technicien d'entretien Thermo Fisher Scientific.

7.21 Préparation du transport du système

AVIS

Les raccordements fluidiques et raccords de capillaires sont extrêmement sensibles à la contamination. La poussière et les débris peuvent contaminer ces raccordements.

Installez toujours des capuchons sur les capillaires et des bouchons sur les raccordements fluidiques ouverts pour éviter toute contamination.

Préparatifs généraux

- 1. Exécutez le script : **B05 Arrêt (de longue durée)** avec le paramètre suivant : **Module à retirer = système entier**.
- Si ce n'est pas déjà fait, verrouillez l'unité d'aiguille (reportez-vous à la section Verrouillage de l'unité d'aiguille (▶ page 251)).
- Pour déplacer légèrement le socle de système et mieux accéder aux pièces du système, déverrouillez le socle du système à l'aide des outils de déverrouillage du socle. Reportez-vous à la section Déverrouillage du socle du système (▶ page 73).
- 4. Placez le système empilé dans une position de travail appropriée.
- 5. Retirez les portoirs d'échantillons et les récipients d'échantillonnage du compartiment à échantillons.

Système de solvant, système de rinçage des joints et système de rinçage de l'aiguille

NOTE Il n'est généralement pas nécessaire de retirer les conduites de solvant des guides de tubulure (même pour remplacer le module à glissière). Toutefois, lorsque vous retirez les conduites de solvant des guides de tubulure dans la pompe et dans tous les modules placés audessus de la pompe dans le système empilé, veillez à ne pas tirer sur les autres tubes présents dans les guides.

1. Videz tous les réservoirs de liquides.

Arrêt

1. Mettez *tous* les modules du système hors tension au moyen de leurs interrupteurs principaux d'alimentation. Il ne suffit pas d'appuyer sur le bouton d'alimentation du système pour couper complètement l'alimentation électrique des modules du système.

NOTE Lorsque le passeur d'échantillon est hors tension, sa porte avant gauche s'ouvre automatiquement afin que le compartiment à échantillons soit correctement ventilé et elle ne peut pas être fermée.

- 2. Débranchez tous les cordons d'alimentation.
- 3. Débranchez tous les câbles de signal vers d'autres appareils.

Nettoyage

- En présence de condensation ou si des échantillons ont été renversés dans le compartiment à échantillons, nettoyez et décontaminez le compartiment à échantillons avant de poursuivre. Vérifiez que le passeur d'échantillon a été nettoyé et/ou décontaminé comme requis.
- 2. Vérifiez si les bacs de fuite contiennent des liquides provenant de fuites. Si c'est le cas, absorbez-les avec une serviette en papier.

Tâches restantes

- 1. Tirez sur les tubes d'évacuation du système pour les faire sortir du bac d'évacuation.
- Fixez les poignées de manutention (reportez-vous à la section Fixation des poignées de manutention (▶ page 350)), le cas échéant.
- 3. Transportez le système.

Redémarrage du système après le transport

- Déverrouillez l'unité d'aiguille (reportez-vous à la section Déverrouillage de l'unité d'aiguille (▶ page 252)).
- Connectez les câbles (reportez-vous à la section Connexion des câbles (▶ page 106)).
- 3. Replacez le système empilé dans la position d'utilisation.
- Verrouillez le socle du système à l'aide des outils de verrouillage (reportez-vous à la section Verrouillage du socle du système (> page 119)).
- 5. Remplissez les réservoirs (reportez-vous aux sections correspondantes).
- Mettez en place les raccordements fluidiques (reportez-vous à la section Mise en place des raccordements fluidiques (> page 120)).
- 7. Mettez le système sous tension (reportez-vous à la section Mise en marche du système (▶ page 148)).

8. Préparez le système pour utilisation (reportez-vous à la section Préparation du système pour l'utilisation initiale (▶ page 139)).

7.22 Fixation des poignées de manutention

À quel moment exécuter cette opération ?

Le système doit être transporté à un autre emplacement.

Éléments nécessaires

4 poignées

Utilisez les poignées de manutention installées lors de l'expédition du système et qui sont munies de 2 tous à vis.



Illustration 228: Poignées de manutention pour l'empilement du système

• 8 vis qui étaient fixées aux poignées

Outils nécessaires

Tournevis type Torx T30

Procédez comme suit



ATTENTION—Lésions corporelles et dommages au système

La fixation des poignées sur le passeur d'échantillon et le levage du système pour causer des lésions corporelles et des dommages au système.

Si vous devez lever et transporter l'ensemble du système, fixez les poignées seulement sur la pompe afin d'éviter de vous blesser et d'endommager le système.

1. Insérez une poignée de manutention dans les rails droite et gauche depuis l'arrière.



Illustration 229: Poignée de manutention glissée sur le rail droit

2. Glissez les poignées vers le devant du système jusqu'à ce que la distance entre le devant du rail et le bord avant de la poignée est d'environ 6 cm.





- 3. Fixez les vis des poignées.
- 4. Insérez une poignée de manutention dans les rails droite et gauche depuis l'arrière.
- 5. Glissez les poignées vers le devant du système pour que le bord arrière des poignées touche le bord arrière des rails.



Illustration 231: Position correcte des poignées de manutention arrière

6. Fixez les vis des poignées.



ATTENTION—Risque de lésions corporelles et de dommages au module

Les poignées risquent de glisser du module si les vis ne sont pas bien fixées. Aussi, le module à glissières peut tomber s'il n'est pas fixé correctement. Cela pourrait causer des lésions corporelles et des dommages au module. Vérifiez les points suivants avant de soulever le système :

- Les poignées sont bien fixées et ne peuvent pas glisser du module.
- Les quatre vis de chaque module à glissières sont bien serrées et aucun module à glissières ne peut tomber (reportez-vous à la section Remplacement du module à glissières (> page 336)).

8 Dépannage

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

8.1 Généralités sur le dépannage

Les fonctions suivantes vous aident à déterminer qu'un problème est survenu pendant l'utilisation du système.

Voyant	Description
Voyants d'état	Le système montre l'état des éléments suivants :
	 Système général : en haut de l'écran dans la Vanquish User Interface. Reportez-vous à la section Icônes d'état de l'état général du système (> page 155).
	 Modules du système Si le microprogramme d'un module du système détecte un problème, un message est affiché dans le volet notification et les voyants d'état suivants du système sont rouges :
	 Barre à LED (reportez-vous à la section Barres à LED (> page 153)) sur le module et représentation de barre à LED dans le volet accueil
	 LED STATUS (reportez-vous à la section Panneaux de commande (> page 149))
Alarmes de fuite	Les fuites constituent un problème potentiel pour la sécurité et la chromatographie. Aussi, si un capteur détecte une fuite, un signal sonore se déclenche afin de vous alerter, en plus du message affiché dans la volet notification, et les voyants d'état deviennent rouges. Suivez les instructions de la section Recherche de fuites de liquide dans le système () page 395).
Messages	Si un microprogramme de module détecte un problème, un message est affiché dans le volet notification. Le volet notification stocke l'information sur tous les avertissements et toutes les erreurs des modules du système et tous les événements du contrôleur du système depuis le dernier démarrage du contrôleur du système.
Problèmes d'analyse	Afin de résoudre les problèmes d'analyse, comme les variations en matière de pics et de temps de rétention, veuillez vous reporter aux sections correspondantes dans Dépannage du système () page 379).

Les sections suivantes contiennent l'information sur les messages et les problèmes de fonctionnement qui peuvent survenir pendant le fonctionnement du système et offrent de l'aide de dépannage.

Aide supplémentaire

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, ayez le numéro de référence et le numéro de série du système à portée de main lorsque vous communiquez avec Thermo Fisher Scientific. Vous trouverez cette information sur la plaque signalétique dans le tiroir du socle du système.

8.2 Messages

Si le microprogramme d'un module du système détecte une erreur, un message est affiché dans le Vanquish User Interface et le système de gestion de données chromatographiques. Les tableaux dans la présente section répertorie les messages les plus courants de chaque module du système et offrent de l'aide 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. Pour limiter la source du problème, du texte supplémentaire peut accompagner un message en particulier, par exemple, dans le système de gestion de données chromatographiques.

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.

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.

8.2.1 Messages communs

La présente section contient de l'information sur les messages communs des modules du système.

Le tableau répertorie les messages les plus courants des modules du système.

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. Si le message s'affiche de nouveau, prenez contact avec l'assistance technique.
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.

Message et code	Description et action corrective
Code 33 Leak detected – eliminate within approx. xx seconds.	Fuite détectée – éliminez la fuite dans les xx secondes suivantes environ.
	xx = le délai maximum en secondes pour résoudre la fuite avant l'arrêt de la séquence
	Trouvez et éliminez la source de la fuite. Si la fuite de la pompe n'est pas réparée dans les 3 minutes, la pompe interrompt le débit.
	 Inspectez visuellement tous les composants et raccords du trajet d'écoulement.
	2. Serrez ou remplacez le raccord ou composant concerné, selon les besoins.
	3. 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.
Code 34 Leak detected.	Fuite détectée.
	Ce message apparaît, car la fuite signalée plus tôt par le code 33 n'a pas été réparée dans les 3 minutes. La pompe s'est arrêtée et/ ou la séquence de fonctionnement a été interrompue à cause de la fuite.
	Trouvez et éliminez la source de la fuite (reportez-vous à la section Recherche de fuites de liquide dans le système () page 395)).
Code 36 Download failed.	Échec du téléchargement.
	Le téléchargement du microprogramme a échoué. Relancez le téléchargement.
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 79 [Device] was not found.	[Appareil] introuvable.
	Le message peut s'afficher pour le capteur de pression dans une tête de pompe ou pour le capteur de pression du système dans le débitmètre (reportez-vous à l'information d'identification dans la section Messages concernant la pompe () page 362)).
	 Tête de pompe : Vérifiez que le câble du capteur de pression de la tête de pompe pour lequel le message s'affiche est correctement branché dans le port de connexion.
	• Débitmètre : Prenez contact avec l'assistance technique.

Message et code	Description et action corrective
Code 84 Could not move X to its destination position. Code 85 Could not move X to its destination position.	 Impossible de mettre X dans la position souhaitée. X = vanne gauche ou droite La vanne de commutation indiquée est peut-être défectueuse. Remplacez-la, reportez-vous à la section Remplacement de la vanne de commutation Vanquish (▶ page 305) ou Remplacement de la vanne de commutation VICI (▶ page 308). 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 – invalid version.	Mauvaise correspondance de microprogramme – version invalide. 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.	 Changement de données pendant la lecture. Tous les modules : Il peut s'agir d'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. En plus, pour la pompe : Si le message s'affiche pour le capteur de pression dans une tête de pompe ou pour le capteur de pression du système dans le débitmètre (reportez-vous à l'information d'identification dans la section Messages concernant la pompe (▶ page 362)) : Tête de pompe : Vérifiez que le câble du capteur de pression de l'appareil indiqué dans le message est correctement branché dans le port de connexion. Débitmètre : Prenez contact avec l'assistance technique

Message et code	Description et action corrective
Code 92 Data verification failed.	Échec de la vérification des données.
	Tous les modules :
	 Il peut s'agir d'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.
	En plus, pour la pompe :
	Si le message s'affiche pour le capteur de pression dans une tête de pompe ou pour le capteur de pression du système dans le débitmètre (reportez-vous à l'information d'identification dans la section Messages concernant la pompe () page 362)) :
	 Tête de pompe : Vérifiez que le câble du capteur de pression de l'appareil indiqué dans le message est correctement branché dans le port de connexion.
	Débitmètre : Prenez contact avec l'assistance technique.
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 traite les données.
	 Dans le système de gestion de données chromatographiques, débranchez et rebranchez le module.
	 Si cela ne résout pas le problème, mettez à jour le firmware ou la version du système de gestion de données chromatographiques.
	3. Si le problème persiste : un logiciel tiers installé sur l'ordinateur, comme les scanners antivirus, ou de mauvaises performances de l'ordinateur peuvent 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.
	2. 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.

Message et code	Description et action corrective
Code 126 Device lost system interlink connection.	L'appareil a perdu la connexion à System Interlink. Tous les autres modules du bus System Interlink sont mis hors tension ou le téléchargement d'un microprogramme est effectué sur l'ensemble des autres modules.
	 Vérifiez si tous les autres modules ont été mis hors tension ou si le téléchargement d'un microprogramme est effectué sur l'ensemble des autres modules du bus.
	 Patientez jusqu'à ce que l'un des modules du bus soit mis sous tension ou jusqu'à ce que le téléchargement du microprogramme soit terminé.
	3 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.
	4. 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.
	2. 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.
Code 131 System interlink bus not	Bus de System Interlink non opérationnel.
operable.	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 133 System Interlink access not	Accès à System Interlink non pris en charge.
supported.	L'accès aux commandes n'est pas permis depuis le contrôleur du système. Vous pouvez uniquement modifier les paramètres du module depuis le système de données chromatographiques.
	Par exemple, attendez la fin de la séquence et que la barre à LED soit verte. Puis, répétez l'opération depuis le contrôleur du système.
Message et code	Description et action corrective
--	--
Code 134 USB access by CDS not	Accès USB par CDS non pris en charge.
supported.	L'accès aux commandes n'est pas permis depuis le système de données chromatographiques. Vous pouvez uniquement modifier les paramètres du module depuis le contrôleur du système.
	Par exemple, attendez la fin du script et que la barre à LED soit verte. Puis, répétez l'opération depuis le système de données chromatographiques.
Code 136 Lock request rejected – already	Requête de verrouillage rejetée – déjà verrouillé par X.
locked by X.	 X = identificateur du support de verrou, avec identificateur des boutons du panneau de commande, adresse USB liée au système de gestion de données chromatographiques 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) ou un bouton du panneau de commande.
	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 identificateur des boutons du panneau de commande, adresse USB liée au système de gestion de données chromatographiques 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 identificateur des boutons du panneau de commande, adresse USB liée au système de gestion de données chromatographiques ou adresse System Interlink liée au contrôleur du système ou à un module
	Le verrou a été libéré automatiquement puisque son support a disparu du System Interlink.
	1. Vérifiez que le module supportant le verrou a été désactivé.
	2. Vérifiez les raccordements System Interlink du module.
Code 152 Assignment or command	Attribution ou commande rejetée – appareil verrouillé.
rejected – device locked.	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).

8.2.2 Messages concernant la pompe

La présente section contient des informations sur les messages avec des numéros de code allant de 4000 à 4999. Ces messages concernent la pompe.

Identification de source supplémentaire

Pour limiter la source du problème, vérifiez le message dans le système de gestion de données chromatographiques pour en savoir plus :

Nom par défaut dans le système de données	Description
PumpModule.Pump	Le message concerne l'ensemble du module de la pompe, y compris le débitmètre (avec le capteur de pression du système).
PumpModule.Pump Left Block (WP)	 Le message concerne les éléments suivants : Bloc gauche de la pompe (A) Composant du bloc gauche de la pompe, par exemple, la tête de pompe Composants du côté gauche de la tête de pompe gauche, par exemple, le capteur de pression à gauche de la tête de pompe ou le piston de travail, livrant %A.
PumpModule.Pump Left Block (EP)	 Le message concerne les éléments suivants : Bloc gauche de la pompe (A) Composant du bloc gauche de la pompe, par exemple, la tête de pompe Composants du côté droit de la tête de pompe gauche, par exemple, le capteur de pression à droite de la tête de pompe ou le piston d'équilibrage, livrant %A
PumpModule.Pump Right Block (WP)	 Le message concerne les éléments suivants : Bloc droite de la pompe (B) Composant du bloc droite de la pompe, par exemple, la tête de pompe Composants du côté gauche de la tête de pompe droite, par exemple, le capteur de pression à gauche de la tête de pompe ou le piston de travail, livrant %B
PumpModule.Pump Right Block (EP)	 Le message concerne les éléments suivants : Bloc droite de la pompe (B) Composant du bloc droite de la pompe, par exemple, la tête de pompe Composants du côté droit de la tête de pompe droite, par exemple, le capteur de pression à droite de la tête de pompe ou le piston d'équilibrage, livrant %B

Messages concernant la pompe

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. Si le message s'affiche de nouveau, prenez contact avec l'assistance technique.
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 4098 Upper pressure limit exceeded	Dépassement de la limite de pression supérieure. Vérifiez si le trajet d'écoulement est obstrué. Exécutez le script : D01 – Tester la contre-pression du système .
Code 4099 Pressure fallen below lower limit.	 Chute de pression en dessous de la limite inférieure. Vérifiez les causes suivantes : Un ou plusieurs réservoirs de solvant sont vides. Exécutez le script : B01 – Modifier les liquides / solvants (Rafraîchir Uniquement = activé). Des bulles d'air sont peut-être piégées dans les conduites de solvant. Vérifiez les filtres de conduite de solvant. Purgez la pompe. Le débit appliqué ne permet pas à la pompe de créer suffisamment de pression pour atteindre la limite de pression inférieure. Modifiez la limite de pression inférieure ou augmentez le débit. Il peut y avoir une fuite dans le système. Exécutez le script : E04 – Test de fuite détaillé (pompe). Un clapet de non-retour peut être contaminé ou défectueux. Nettoyez ou remplacez les clapets de non-retour. Reportez-vous à la section Entretien des clapets de non-retour de la tête de pompe () page 226).
Code 4100 Purge pressure limit exceeded	 Dépassement de la limite de pression de purge Génération d'une pression supérieure à 5 MPa (50 bar) après lancement d'un cycle de purge. Exécutez le script : D01 – Tester la contre-pression du système. Prenez les mesures correctives suggérées par le script. Si le message s'affiche avec un message concernant la contre- pression du filtre du débitmètre (code 4533), remplacez le filtre (reportez-vous à la section Remplacement du filtre du débitmètre (▶ page 232)). Une vanne de purge peut être défectueuse. Prenez contact avec l'assistance technique.

Le tableau répertorie les messages les plus courants de la pompe.

Message et code	Description et action corrective
Code 4101 Not ready. Please run self- test.	Non prêt. Veuillez exécuter un test automatique. La pompe n'est pas prête, car le test automatique a échoué. Répétez le test automatique. Exécutez le script : B04 – Test automatique du système .
Code 4112 Self-test failed because of fluctuating pressure.	Échec du test automatique en raison de variations de pression. Une fluctuation de pression s'est produite lors du test automatique. Patientez jusqu'à ce que la pression du système soit redescendue à zéro, ou dépressurisez le système et répétez le test automatique. Exécutez le script : B04 – Test automatique du système .
Code 4113 Self-test failed because pressure is out of range.	Échec du test automatique en raison d'une pression hors plage. La pression n'est pas redescendue à zéro lors du test automatique. Vérifiez que les câbles de capteur de pression sont correctement branchés. Patientez jusqu'à ce que la pression du système soit redescendue à zéro, ou dépressurisez le système et répétez le test automatique. Exécutez le script : B04 – Test automatique du système . Si le message s'affiche de nouveau, le capteur de pression du système est peut-être défectueux. Prenez contact avec l'assistance technique.
Code 4114 The device is busy. Please retry after 1 minute.	L'appareil est occupé. Veuillez essayer dans 1 minute. La pompe n'est pas prête, par exemple en raison d'un test automatique ou d'une procédure de ré-étalonnage en cours. Réessayez dans 1 minute. Si le message s'affiche de nouveau, patientez jusqu'à ce que la pression du système s'abaisse ou dépressurisez le système, puis éteignez et rallumez la pompe.
Code 4119 Can't start pump while alarm is on.	Impossible de démarrer la pompe tant qu'une alarme est en cours. Une alarme est en cours, par exemple suite à la détection d'une fuite. Vous pouvez relancer le débit de pompe uniquement lorsque l'alarme a été résolue. Désactivez le signal sonore de l'alarme en appuyant sur le bouton MUTE ALARM sur le panneau de commande de la pompe. Trouvez et éliminez la source de l'alarme en suivant les instructions fournies pour le message qui accompagne l'alarme.

Message et code	Description et action corrective
Code 4120 The rear seal wash system has run out of wash solution.	Le liquide de rinçage du système de rinçage des joints arrière est épuisé.
	Ce message vous avertit uniquement si la fonction de détecteur de gouttes du détecteur de rinçage des joints est activée (réglage standard).
	 Le réservoir de liquide de rinçage des joints peut être vide. Ajoutez du liquide frais de rinçage des joints et exécutez le script : B01 – Modifier les liquides / solvants (Liquide de rinçage des joints arrière, Rafraîchir Uniquement = activé). Respectez les recommandations concernant le liquide de rinçage de des joints de la section Remplissage du réservoir de liquide de rinçage des joints (▶ page 134).
	 Vérifiez si les conduites de rinçage des joints sur l'ensemble du trajet d'écoulement présentent des signes de fuite ou d'obstruction. Toutes les conduites doivent être correctement connectées et acheminées. Remplacez les conduites de rinçage des joints selon les besoins. Reportez-vous à la section Entretien du système de rinçage des joints (> page 282).
	 Vérifiez si du liquide s'échappe des raccordements du système de rinçage des joints ou sous les têtes de pompe. Reportez-vous à la section Test d'étanchéité du système de rinçage des joints (> page 282).
	 Vérifiez si les tubes péristaltiques présentent des signes d'obstruction ou de détérioration. Remplacez les tubes selon les besoins. Reportez-vous à la section Remplacement du tube péristaltique dans la pompe de rinçage des joints (> page 293).
	 Vérifiez que le tube péristaltique est correctement inséré dans la pompe de rinçage des joints et que le levier de la pompe n'est pas bloqué.
	 Vérifiez si le tube de rinçage des joints est connecté au bon orifice dans le détecteur de rinçage des joints (orifice étiqueté in ; reportez-vous à l'image dans la section Test d'étanchéité du système de rinçage des joints (> page 282)).
	 Vérifiez si les électrodes du détecteur de rinçage des joints sont contaminées ou détériorées. Remplacez le détecteur de rinçage des joints selon les besoins. Reportez-vous à la section Remplacement du détecteur de rinçage des joints (> page 298).
Code 4121 Piston seal leakage has	La fuite du joint de piston dépasse la limite recommandée.
exceeded the recommended limit.	Ce message vous avertit uniquement si la fonction de détecteur de gouttes du détecteur de rinçage des joints est activée (réglage standard).
	Vérifiez s'il y a des fuites dans la pompe. Exécutez le script : E04 – Test de fuite détaillé (pompe) . Quand les joints de piston fuient, remplacez-les. Reportez-vous à la section Remplacement du joint du piston () page 209).

Message et code	Description et action corrective
Code 4122 The rear seal leak sensor detects drops constantly.	Le capteur de fuite de joint arrière détecte constamment des gouttes.
	Ce message vous avertit uniquement si la fonction de détecteur de gouttes du détecteur de rinçage des joints est activée (réglage standard).
	 Vérifiez que le liquide s'écoule correctement dans le dispositif d'évacuation par l'orifice d'évacuation situé en bas à droite sur le bac de fuite.
	 Inspectez la face inférieure du détecteur de rinçage des joints : Retirez le bloc de détecteur de rinçage des joints de la pompe et déconnectez le détecteur du pied. Séchez la face inférieure avec un chiffon ou une serviette en papier. Inspectez les électrodes du détecteur. Si elles sont propres et en bon état, ré-assemblez le détecteur et le pied et réinstallez le bloc dans la pompe. Reportez-vous à la section Remplacement du détecteur de rinçage des joints (▶ page 298).
	 Si le message s'affiche de nouveau, le détecteur de rinçage des joints peut être défectueux. Remplacez le détecteur de rinçage des joints. Reportez-vous à la section Remplacement du détecteur de rinçage des joints (> page 298).
Code 4125 Degasser malfunction	Dysfonctionnement du dégazeur
	La fonction de surveillance du niveau de vide détecte un vide insuffisant au niveau du dégazeur.
	Mettez à jour le micrologiciel vers la version 2.41 ou ultérieure.
Code 4127 The pump drive is still in undock position. Please execute a dock command.	La commande de la pompe est toujours en position de retrait. Veuillez exécuter une commande Dock.
	Ce message vous avertit lors la maintenance de la pompe (par exemple, lors du remplacement de pistons ou des joints de piston) si vous essayez de démarrer la pompe alors que la tête de pompe n'est pas encore dans la position appropriée pour une utilisation normale.
	Replacez la tête de pompe dans la position d'utilisation : sur le panneau de commande, les LED situées à côté du bouton FLOW indiquent quelle tête de pompe est dans la position de maintenance (la LED clignote en vert). Vérifiez que cette tête de pompe est sélectionnée (la LED située à côté du bouton SELECT est verte) et appuyez sur le bouton DOCK . Lorsque la pompe est en position d'utilisation, la LED située à côté du bouton FLOW arrête de clignoter.
Code 4148 Can't perform this command while the flow is on.	Impossible d'exécuter cette commande alors que le débit est activé. Vous avez lancé un test automatique alors que le débit n'était pas
	redescendu à zéro. Arrêtez le débit et répétez la commande.

Message et code	Description et action corrective
Code 4152 Pressure sensor malfunction. Check cable connection and retry.	Dysfonctionnement du capteur de pression. Vérifiez la connexion de câble et réessayez.
	Vérifiez si le câble du capteur de pression pour lequel le message s'affiche est correctement branché dans le port de connexion. Redémarrez le débit.
	Si le message s'affiche de nouveau, le capteur de pression peut être défectueux :
	 Si le message apparaît pour le capteur de pression d'une tête de pompe, remplacez la tête de pompe.
	 Si le message apparaît pour le capteur de pression de système, prenez contact avec l'assistance technique.
Code 4158 Out of eluent X.	Éluant X épuisé.
	Ce message vous avertit uniquement si vous surveillez la consommation de solvant. Il indique le solvant dont le réservoir est vide. Remplissez le solvant. Exécutez le script : B01 – Modifier les liquides / solvants .
Code 4161 Pressure recalibration	Le réétalonnage de la pression diverge de xx bar.
deviates by xx bar.	La pression de la pompe n'était pas redescendue à zéro lors du test automatique. Patientez jusqu'à ce que la pression soit redescendue, puis répétez le test. Exécutez le script : B04 – Test automatique du système .
Code 4208 System pressure too high. Please relieve pressure and retry.	Pression du système trop élevée. Veuillez relâcher la pression et réessayer.
	Procédez comme suit :
	 Vérifiez que les câbles de capteur de pression sont correctement raccordés. Si nécessaire, rebranchez les câbles.
	 Vérifiez si le trajet d'écoulement est obstrué. Exécutez le script : D01 – Tester la contre-pression du système.
	 Effectuez un test automatique. Exécutez le script : B04 – Test automatique du système.
	Si le message s'affiche de nouveau, le capteur de pression du système est peut-être défectueux. Prenez contact avec l'assistance technique.
Code 4209 Could not build up enough pressure. Please check the fitting plug.	Impossible de générer une pression suffisante. Veuillez vérifier le bouchon de raccordement.
	La pompe n'a pas pu générer assez de pression lors de l'étalonnage du capteur de pression. Cela peut être dû à :
	 De l'air potentiellement emprisonné dans le système. Purgez la pompe. Exécutez le script : C02 – Purger la pompe.
	 Il est possible qu'une fuite soit présente. Exécutez le script : E04 Test de fuite détaillé (pompe).

Message et code	Description et action corrective
Code 4213 Head pressure does not match system pressure. Execute the self- test or perform pressure recalibration.	La pression de tête ne correspond pas au signal de pression détecté par le système. Exécutez le test automatique ou effectuez un ré-étalonnage de la pression.
	 Vérifiez que le trajet d'écoulement depuis la tête de la pompe à l'orifice d'admission du débitmètre ne présente aucun signe de blocage, en particulier les extrémités Viper. Si vous suspectez un blocage, remplacez le capillaire.
	 Effectuez un test automatique. Exécutez le script : B04 – Test automatique du système.
	 Si le message persiste, effectuez un nouvel étalonnage de la pression. Exécutez le script : C05 – Régler les capteurs de pression du système.
	4. Si le message s'affiche de nouveau, prenez contact avec l'assistance technique.
Code 4220 Pump head not recognized. Please plug in the pressure sensor cable.	Tête de pompe non reconnue. Veuillez brancher le câble du capteur.
	 Branchez le câble du capteur de pression de la tête de pompe sur le connecteur Pressure. Si le câble est branché, vérifiez s'il est bien inséré.
	2. Si le problème persiste, la tête de pompe est peut-être défectueuse. Prenez contact avec l'assistance technique ou remplacez la tête de pompe (reportez-vous à la section Remplacement de la tête de pompe (▶ page 197)).
Code 4231 Pressure sensor not recognized. Please plug in the sensor cable.	Capteur de pression non reconnu. Veuillez brancher le câble du capteur.
	Vérifiez si le câble du capteur de pression pour lequel le message s'affiche est correctement branché dans le port de connexion. Redémarrez le débit.
	Si le message s'affiche de nouveau, le capteur de pression peut être défectueux :
	 Si le message apparaît pour le capteur de pression d'une tête de pompe, remplacez la tête de pompe (reportez-vous à la section Remplacement de la tête de pompe (> page 197)).
	 Si le message s'affiche pour le capteur de pression du système, il est peut-être défectueux. Prenez contact avec l'assistance technique.
Code 4232 The identification number of	Le numéro d'identification de XX a été modifié.
the XX has changed.	XX est la tête de pompe gauche ou la tête de pompe droite
	Ce message s'affiche lorsqu'une tête de pompe a été remplacée ou lorsque les têtes de pompe gauche et droite ont été interverties, par exemple de façon accidentelle pendant une opération de maintenance ou d'entretien. Ce message est affiché à titre d'information uniquement. Aucune action n'est requise.
Code 4234 The flow was stopped due to a request from another module.	Le débit a été interrompu suite à une requête provenant d'un autre module.
	Le débit de la pompe a été arrêté en raison d'un problème avec un autre module. Vérifiez les messages des autres modules.

Message et code	Description et action corrective
Code 4238 Mainboard initialization failed.	Échec de l'initialisation de la carte mère. L'initialisation de la carte mère a échoué. 1. Cela peut être une erreur temporaire. Mettez le module hors tension. Patientez 5 secondes, puis remettez le module sous tension. 2. Si le message s'affiche à nouveau, mettez à jour le microprogramme. 3. Si le problème persiste, la carte mère est peut-être défectueuse. Prenez contact avec l'assistance technique.
Code 4503 The pressure exceeded the absolute limit.	La pression dépasse la limite absolue. Vérifiez si le trajet d'écoulement est obstrué. Exécutez le script : D01 – Tester la contre-pression du système .
Code 4504 NC drive initialization error.	Erreur d'initialisation de l'entraînement NC. Répétez le test automatique. Exécutez le script : B04 – Test automatique du système.
Code 4508 Piston drive home sensor remains dark.	Le capteur d'entraînement du piston reste éteint. Répétez le test automatique. Exécutez le script : B04 – Test automatique du système .
Code 4512 Cannot pressurize due to inlet valve failure or air in system.	 Impossible de pressuriser en raison d'une défaillance de la vanne d'entrée ou de la présence d'air dans le système. Il a fallu répéter une phase de compression, puisque la pression attendue n'a pas été atteinte. Cela peut être causé par une présence d'air dans le système ou parce que le clapet de nonretour d'aspiration doit être nettoyé. Toutefois, comme la phase de compression suivante était réussie, la pompe livre le bon débit. Si le message est rare et qu'il n'est pas accompagné d'autres messages, par exemple code 4532 ou code 4538, vous pouvez le tolérer. Autrement : Contrôlez le niveau de liquide dans le réservoir de solvant. Purgez la pompe. Exécutez le script : C02 – Purger la pompe. Si le message s'affiche de nouveau après la purge, nettoyez le clapet de non-retour d'aspiration (reportez-vous à la section Nettoyage des clapets de non-retour (tête de pompe) (▶ page 229)). Utilisez seulement les véritables conduites de solvant autorisées par Thermo Fisher Scientific pour cette pompe.
Code 4513 Home position deviation XX. Please execute self-test.	Déviation de la position initiale XX. Veuillez exécuter un test automatique. XX est la valeur de la déviation. Assurez-vous d'utiliser le liquide de rinçage des joints recommandé (reportez-vous à la section Remplissage du réservoir de liquide de rinçage des joints () page 134)) et effectuez un test automatique. Exécutez le script : B04 – Test automatique du système.

Message et code	Description et action corrective
Code 4514 The user defined calibration data of the flow sensor is invalid.	Les données d'étalonnage du capteur de débit ne sont pas valides. Les données d'étalonnage du solvant sélectionné sont incomplètes. Par exemple, parce que la procédure d'étalonnage a été interrompue de manière inattendue. Répétez l'étalonnage du solvant sélectionné (reportez-vous à la section Ajout de solvants personnalisés au tableau d'étalonnage (▶ page 170)).
Code 4517 Residual pressure too high for flow autozero.	Pression résiduelle trop élevée pour remise automatique du débit à zéro. Arrêtez le débit et patientez jusqu'à ce que la pression soit descendue.
Code 4519 Calibration table full.	Tableau de calibration plein. Supprimez un solvant du tableau d'étalonnage de solvants pour libérer de l'espace. Dans le volet script , allez à Étalonnage du solvant . Exécutez le script : M34 – Supprimer les données d'étalonnage du solvant.
Code 4520 Flowmeter memory unavailable.	Mémoire du débitmètre indisponible. Vérifiez si le débitmètre est correctement inséré.
Code 4522 Unknown solvent.	Solvant inconnu. Le solvant sélectionné n'est pas disponible dans le tableau d'étalonnage des solvants. Cela peut être dû aux raisons suivantes : Le solvant à été retiré du tableau de calibration. Le débitmètre a été remplacé et un solvant (du même nom) n'a pas été étalonné et ajouté à la configuration du nouveau débitmètre. Sélectionnez un solvant dans le tableau d'étalonnage de solvants ou ajouter le tableau d'étalonnage des solvants (reportez-vous à la section Ajout de solvants personnalisés au tableau d'étalonnage (▶ page 170)).
Code 4523 Drift rate too high for flow autozero.	 Taux de dérive trop élevé pour remise automatique du débit à zéro. Le débit n'est pas stable. 1. Patientez jusqu'à ce que la pression du système redescende à zéro. 2. Purgez la pompe pour éliminer tout air résiduel. 3. Quand la pompe est sous tension depuis au moins 6 heures, répétez la remise automatique à zéro.
Code 4526 Command sequence error.	Erreur de séquence de commandes. Les commandes ont été exécutées dans un ordre erroné ou inattendu. Recherchez des erreurs dans votre méthode de script de commandes.
Code 4529 The pump capillary connections are setup for solvent calibration. Normal operation is not possible.	Les connexions capillaires de la pompe sont configurées pour l'étalonnage du solvant. Le fonctionnement normal n'est pas possible. Une tentative d'étalonnage a été interrompue. Utilisez les scripts d'étalonnage de solvants pour complétez l'étalonnage ou retournez la pompe à la configuration capillaire d'origine et raccordez tous les capillaires et le système selon les besoins de l'application.

Message et code	Description et action corrective
Code 4531 Cannot deliver flow at this pressure.	 Impossible de livrer un débit à cette pression. La pompe ne peut pas livrer le débit demandé à la pression résultante. Adaptez la configuration de l'application. Par exemple, diminuez le débit. Reportez-vous à la section Interdépendance entre la pression de la pompe et le débit (> page 169).
Code 4532 Cannot regulate flow due to air in system or inlet valve failure.	 Impossible de réguler le débit en raison d'une présence d'air dans le système ou d'un défaut de la vanne d'admission. Une ou plusieurs phases de compression n'ont pas abouti. La pompe n'a pas pu livrer le début adéquat. Cela peut être causé par une présence d'air dans le système ou parce que le clapet de non-retour d'aspiration doit être nettoyé. Le message apparaît avec le message code 4512. Contrôlez le niveau de liquide dans le réservoir de solvant. Purgez la pompe. Exécutez le script : CO2 – Purger la pompe. Si le message s'affiche de nouveau après la purge, nettoyez le clapet de non-retour d'aspiration (reportez-vous à la section Nettoyage des clapets de non-retour (tête de pompe) () page 229)).
Code 4533 Elevated filter backpressure. Consider replacing the flowmeter filter soon.	 Utilisez seulement les véritables conduites de solvant autorisées par Thermo Fisher Scientific pour cette pompe. Contrepression élevée dans le filtre. Envisagez de remplacer bientôt le filtre du débitmètre. Ce message apparaît pour vous signaler un problème en cours. Le message indique qu'il peut y avoir un problème avec les solvants, par exemple l'apparition d'algues. Envisagez de remplacer le filtre du débitmètre (reportez-vous à la section Remplacement du filtre du débitmètre (> page 232)). Respectez les recommandations de la section Utilisation de solvants et d'additifs (> page 145).
Code 4534 System pressure exceeded maximum pressure during purge.	 La pression du système dépasse la pression maximale pendant une purge. 1. Sur la sortie du débitmètre, retirez le capillaire vers le passeur d'échantillon et réessayez. 2. Dépend des éléments suivants : L'erreur n'est plus présente : Les fluidiques du passeur d'échantillon ou le capillaire vers le passeur d'échantillon peuvent être obstrués. Selon les besoins, éliminez la source de l'obstruction et/ou remplacez le capillaire entre le débitmètre et le passeur d'échantillon. L'erreur est toujours présente : dans le débitmètre, l'élément en Y ou le capillaire de restriction sont peut-être obstrués. Prenez contact avec l'assistance technique.
Code 4536 Solvent calibration failed.	Échec de l'étalonnage du solvant. Raccords de capillaires ou bouchons de la tête de pompe incorrectement installés ou insuffisamment serrés. 1. Si nécessaire, reconnectez ou resserrez. 2. Renouvelez l'étalonnage.

Message et code	Description et action corrective
Code 4538 Flow too low – check for leaks.	Débit trop faible – vérifiez l'absence de fuites.
	apparaît seul ou avec le code de message 4512 .
	• Le message est accompagné du code de message 4512 : effectuez les actions correctives suggérées pour le message d'accompagnement. Si cela ne règle pas le problème, il peut y avoir une fuite dans le système. Trouvez et éliminez la source de la fuite (reportez-vous à la section Recherche de fuites de liquide dans le système () page 395)).
	 Le message apparaît seul : Il peut y avoir une fuite dans le système. Trouvez et éliminez la source de la fuite (reportez-vous à la section Recherche de fuites de liquide dans le système (> page 395)).
Code 4539 Flow too high.	Débit trop élevé.
	La pompe ne peut pas réduire le débit au débit nominal.
	 Des messages apparaissent lorsque le débit configuré est proche de zéro (par exemple, lors de l'arrêt du débit de la pompe) : ajustez les capteurs de pression et de débit. Exécutez les scripts : C05 – Régler les capteurs de pression du système et C01 – Régler les décalages du capteur de débit de la pompe.
	 Un message apparaît dans d'autres situations : contactez l'assistance technique.

8.2.3 Messages concernant le passeur d'échantillon

La présente section contient de l'information sur les messages avec des numéros de code entre 5000 et 5999. Ces messages concernent le passeur d'échantillon.

Le tableau répertorie les messages les plus courants du passeur d'échantillon.

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. Si le message s'affiche de nouveau, prenez contact avec l'assistance technique.
Code 5002 Command rejected – module is in error state.	Commande rejetée - le module est en état d'erreur. Une erreur interne est survenue. Effectuez un test automatique.
Code 5010 Horizontal needle drive blocked. Loosen transport lock completely.	Entraînement horizontal de l'aiguille bloqué. Desserrez complètement le verrouillage d'expédition. Desserrez la vis de verrouillage d'expédition au-dessus de l'unité d'aiguille. Reportez-vous à la section Déverrouillage de l'unité d'aiguille (> page 252). Vérifiez si l'entraînement horizontal de l'aiguille peut être déplacé aisément. Retirez tout obstacle présent sur le trajet de l'entraînement de l'aiguille.
Code 5011 Vertical needle drive blocked. Check for correct installation of sample loop.	Entraînement vertical de l'aiguille bloqué. Vérifiez l'installation de la boucle d'échantillonnage. Vérifiez si la boucle d'échantillonnage est correctement installée. Reportez-vous aux étapes d'installation de la boucle d'échantillonnage de la section.
Code 5012 Selftest failed.	Échec du test automatique. Le test automatique du passeur d'échantillon a échoué. Vérifiez s'il y a des messages supplémentaires indiquant quel est le problème. Prenez contact avec l'assistance technique de Thermo Fisher Scientific pour obtenir de l'aide.
Code 5013 Carousel blocked. Check for obstructions with racks or sample containers.	Carrousel bloqué. Recherchez d'éventuelles obstructions par des portoirs d'échantillons ou des récipients d'échantillonnage. Faites pivoter le carrousel et vérifiez s'il se déplace aisément. Retirez tout obstacle présent sur le trajet du carrousel.
Code 5014 Rack not found. Check for correct rack position in the carousel and sample sequence.	Portoir introuvable. Vérifiez que le portoir est correctement positionné dans le carrousel et la séquence de prélèvement des échantillons. Vérifiez si un portoir d'échantillons ou une plaque à puits est installé dans ce segment du carrousel. Vérifiez s'il est placé correctement (reportez-vous à la section) et corrigez sa position si nécessaire.

Message et code	Description et action corrective
Code 5015 Vial not found. Check for correct vial position in the rack and sample sequence.	Échantillon introuvable. Vérifiez que l'échantillon est correctement positionné dans le portoir et la séquence de prélèvement des échantillons.
	Vérifiez si un flacon ou une plaque à puits est présent à la position indiquée. Placez un récipient d'échantillonnage à la position spécifiée (reportez-vous à la section). Vérifiez la position des échantillons définie dans la séquence de prélèvement des échantillons. Spécifiez une position différente, si nécessaire.
Code 5016 Needle seat not found. Install needle seat.	Siège d'aiguille introuvable. Installez le siège d'aiguille. Vérifiez si le siège d'aiguille est installé. Installez ou remplacez le siège d'aiguille et l'unité d'aiguille, selon les besoins (reportez- vous à la section Remplacement du siège de l'aiguille et de l'unité de l'aiguille () page 254)).
Code 5017 Needle hit an unexpected	L'aiguille a heurté un objet inattendu en raison d'une configuration
well plate configuration.	Vérifiez si la configuration de portoir d'échantillons ou de la plaque à puits puits est correctement définie et si la position des échantillons est appropriée dans la séquence. Prenez en considération le nombre de positions du portoir d'échantillons ou de la plaque à puits configuré, ainsi que la hauteur de plaque à puits correspondante.
Code 5018 Vial pusher hit an object that is higher than expected. Check the rack and/or well plate configuration.	Le pousseur de flacon a heurté un objet de hauteur plus importante que prévu. Vérifiez la configuration du portoir d'échantillons et / ou de la plaque à puits.
	Assurez-vous que le bon portoir d'échantillons ou la bonne plaque à puits sont configurés. Prenez en considération la hauteur du portoir d'échantillons ou de la plaque à puits.
Code 5020 Septum too strong/tight for bottom detection. Check documentation for supported consumables.	Septum trop important / serré pour une détection du fond. Consultez la documentation pour connaître les consommables pris en charge.
	La détection du fond n'est pas opérationnelle avec le septum actuel. Utilisez un septum pris en charge ou désactivez la détection du fond. Pour les capuchons d'échantillons et les couvercles de plaque à puits, reportez-vous à la section Recommandations sur les bouchons de flacons et les couvercles des plaques à puits () page 167).
Code 5021 Insulation cover removed while temperature control is on. Replace cover or temperature control will shut down in X.X minutes.	Couvercle isolant retiré pendant la régulation de la température. Remettez en place le capot ou la régulation de la température s'arrêtera dans X.X minutes.
	Où X.X = le délai en minutes
	Pour les passeurs d'échantillon munis d'un couvercle isolant (boucle), celui-ci doit être installé pour utiliser la thermostatisation du compartiment à échantillons. En cas d'installation inappropriée, la thermostatisation ne pourra être activée ou sera arrêtée automatiquement afin de protéger les composants internes du taux d'humidité élevé et de la forte condensation qui peut apparaître dans le compartiment à échantillons lors de la thermostatisation.
	Si le message apparaît alors que le couvercle est monté, vérifiez que la partie supérieure du couvercle (où se trouve l'aimant) est solidement installée. Vous pouvez surveiller dans l'Instrument Audit Trail si le couvercle est correctement détecté lorsque vous le retirez et l'installez.

Message et code	Description et action corrective
Code 5022 Temperature control shut down due to missing loop cover.	Arrêt de la régulation de la température en raison de l'absence du capot de boucle. Pour les passeurs d'échantillon munis d'un couvercle isolant (boucle), celui-ci doit être installé pour utiliser la thermostatisation du compartiment à échantillons. En cas d'installation inappropriée, la thermostatisation ne pourra être activée ou sera arrêtée automatiquement afin de protéger les composants internes du taux d'humidité élevé et de la forte condensation qui peut apparaître dans le compartiment à échantillons lors de la thermostatisation. Si le message apparaît alors que le couvercle est monté, vérifiez que la partie supérieure du couvercle (où se trouve l'aimant) est solidement installée. Vous pouvez surveiller dans l'Instrument Audit Trail si le couvercle est correctement détecté lorsque vous le retirez et l'installez.
Code 5023 Insulation Loop Cover Bracket removed while temperature control is on. Replace frame or temperature control will shut down in X.X minutes.	Retrait du support du couvercle isolant de la boucle pendant le fonctionnement du contrôle de la température. Remettez-le en place, sinon la régulation de la température s'arrêtera dans X.X minutes. Où X.X = le délai en minutes Le support de montage du couvercle isolant a été retiré alors que la régulation de la température était activée. Pour utiliser la fonctionnalité de thermostatisation du compartiment à échantillons, le couvercle isolant doit être installé. Réinstallez ou remplacez le support de montage du couvercle isolant comme décrit à la section Remplacement du siège de l'aiguille et de l'unité de l'aiguille () page 254) et installez le couvercle isolant comme décrit à la section Installation du couvercle isolant () page 98).
Code 5024 Insulation Loop Cover Bracket missing.	Support du couvercle isolant de la boucle manquant. Le support de montage du couvercle isolant a été retiré alors que le contrôle de la température était activé. Réinstallez ou remplacez le support de montage du couvercle isolant comme décrit à la section Remplacement du siège de l'aiguille et de l'unité de l'aiguille (▶ page 254) et installez le couvercle isolant comme décrit à la section Installation du couvercle isolant (▶ page 98).
Code 5029 Carousel home sensor failure.	Échec capteur du carrousel. Lorsque le message s'affiche, notez son code, puis mettez le module hors tension. Patientez 5 secondes, puis remettez le module sous tension. Si le message s'affiche de nouveau, prenez contact avec l'assistance technique.
Code 5030 Carousel motor failure or blockade.	Échec capteur du carrousel. Lorsque le message s'affiche, notez son code, puis mettez le module hors tension. Patientez 5 secondes, puis remettez le module sous tension. Si le message s'affiche de nouveau, prenez contact avec l'assistance technique.

Message et code	Description et action corrective
Code 5032 Both doors not closed while temperature control is on. Close both doors or temperature control will shut	Portes non fermées alors que la régulation de la température est activée. Fermez les portes, sinon la régulation de la température s'arrêtera dans xx minutes.
down in xx minutes.	xx = le nombre de secondes avant l'arrêt du passeur d'échantillon
	Gardez les deux portes du passeur d'échantillon fermées quand la régulation de la température est activée.
Code 5033 Temperature control shut down due to open door.	Arrêt de la régulation de la température en raison de l'ouverture de la porte.
	Fermez les deux portes du passeur d'échantillon et réactivez la régulation de la température.
Code 5035 Temperature control shut down due to xx fan failure.	Arrêt de la régulation de la température en raison d'une défaillance de ventilation xx.
	xx = le ventilateur touché
	Lorsque le message s'affiche, notez son code, puis mettez le module hors tension. Patientez 5 secondes, puis remettez le module sous tension. Si le message s'affiche de nouveau, prenez contact avec l'assistance technique.
Code 5037 Backup memory error.	Erreur mémoire de sauvegarde. Module non opérationnel.
Module not operational.	Lorsque le message s'affiche, notez son code, puis mettez le module hors tension. Patientez 5 secondes, puis remettez le module sous tension. Si le message s'affiche de nouveau, prenez contact avec l'assistance technique.
Code 5039 Calibration process refused,	Processus d'étalonnage refusé, pression en dessous de 100 bars.
pressure below 100 bar.	Il est impossible d'effectuer l'étalonnage du capteur de pression, car la pression du système est trop basse.
	 Inspectez visuellement le système pour voir s'il présente des fuites, si des pièces sont endommagées ou si certains raccordements sont ouverts.
	 Purgez le passeur d'échantillon : exécutez le script : C04 – Purger l'échantillonneur.
	 Vérifiez s'il y a des fuites ou de l'air dans le système : exécutez le script : D02 – Tester le serrage du système (ce qui devrait être testé = l'échantillonneur).
	 Répétez l'étalonnage du capteur de pression : exécutez le script : C05 – Régler les capteurs de pression du système.
Code 5040 Model – Firmware Mismatch! This firmware only supports VN-A10 models! Download appropriate	Modèle – Microprogramme inadéquat ! Ce microprogramme ne prend en charge que les modèles VN-A10 ! Téléchargez le microprogramme approprié !
firmware!	Vous avez essayé de télécharger le mauvais microprogramme. Sélectionnez le fichier du microprogramme du passeur d'échantillon Vanquish Neo pour le téléchargement.
Code 5041 Abort error occurred.	Erreur d'abandon. Séquence de commandes abandonnée.
Command sequence aborted.	Vérifiez s'il y a des messages supplémentaires indiquant quel est le problème.

Message et code	Description et action corrective
Code 5042 Vanquish Neo sampler detected unknown inject valve bypass position. Set correct position and run self-test command.	L'échantillonneur Vanquish Neo a détecté une position de bypass de vanne d'injection inconnue. Réglez-la à la position adéquate et exécutez la commande de test automatique. Le passeur d'échantillon a été interrompu pendant un script, comme A06 – Modifier le circuit fluidique / flux de tâches . Répétez le script.
Code 5043 Vanquish Neo Pump not found on system interlink. Cannot receive pressure from pump SN xx.	Pompe Vanquish Neo introuvable sur System Interlink. Impossible de recevoir la pression depuis la pompe NS xx. Où xx = le numéro de série de la pompe Vérifiez si tous les câbles System Interlink du système Vanquish sont correctement branchés.
Code 5044 Validation of vial bottom detection failed.	Échec de validation de la détection du fond de flacon. Exécutez la commande SelfTest (de test automatique) ou mettez hors tension le passeur d'échantillon, puis mettez-le à nouveau sous tension pour répéter la validation.
Code 5045 Could not pressurize while delivering solvent. Check for air or leakages in system.	Pressurisation impossible pendant la distribution du solvant. Vérifiez qu'il n'y a pas d'air ou de fuites dans le système. Exécutez le script suivant et suivez les instructions : D02 – Tester le serrage du système (ce qui devrait être testé = l'échantillonneur) .
Code 5099 The device is busy.	L'appareil est occupé. Le passeur d'échantillon n'est pas prêt, par exemple parce qu'un test automatique est en cours ou parce qu'il est toujours en train d'exécuter une autre commande. Essayez à nouveau lorsque le passeur d'échantillon est prêt.

8.2.4 Messages concernant le compartiment à colonnes

La présente section contient de l'information sur les messages avec des numéros de code entre 2000 et 2999. Ces messages concernent le compartiment à colonnes.

Le tableau répertorie les messages les plus courants du compartiment à colonnes.

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. Si le message s'affiche de nouveau, prenez contact avec l'assistance technique.
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 2095 Wrong right valve pod installed.	Mauvaise capsule de vanne droite installée. Vérifiez si la capsule de vanne configurée et la capsule de vanne installée correspondent.
Code 2096 Wrong left valve pod installed.	Mauvaise capsule de vanne gauche installée. Vérifiez si la capsule de vanne configurée et la capsule de vanne installée correspondent.
Code 2097 Right valve position error.	 Erreur de position de la vanne droite. 1. Vérifiez si la capsule de vanne configurée et la capsule de vanne installée correspondent. 2. Vérifiez si la capsule de vanne est correctement installée. 3. Lancez la fonction de vanne de captage en appuyant sur le bouton VALVE RIGHT sur le panneau de commande et en le maintenan enfoncé.
Code 2098 Left valve position error.	 Erreur de position de la vanne gauche. 1. Vérifiez si la capsule de vanne configurée et la capsule de vanne installée correspondent. 2. Vérifiez si la capsule de vanne est correctement installée. 3. Lancez la fonction de vanne de captage en appuyant sur le bouton VALVE LEFT sur le panneau de commande.

8.3 Dépannage du système

La présente section recense les conseils à suivre en cas d'incidents pendant l'utilisation du système Vanquish Neo.

Elle fournit des renseignements sur les symptômes et leurs causes concernant directement les modules du système LC et les connexions entre les modules. Pour obtenir de l'information sur la chromatographie en général et le dépannage des applications, consultez la documentation technique.

La Vanquish User Interface offre quelques scripts de dépannage du système.

Problèmes de détection

Problème	Script
 Contre-pression élevée dans la colonne Arrêt pour cause de surpression Mauvaise précision de l'aire des pics Augmentation continuelle de la pression Aucun débit ou la pompe a émis un message indiquant qu'elle ne pouvait pas réguler le débit 	D01 – Tester la contre-pression du système
 Faible pression dans la colonne Variations dans les temps de rétention Mauvaise précision de l'aire des pics Aucun débit ou la pompe a émis un message indiquant qu'elle ne pouvait pas réguler le débit 	D02 – Tester le serrage du système (ce qui devrait être testé = le système)

Problèmes pendant l'analyse

Problème	Script
L'injection ou la séquence se sont arrêtées et le script B03 – Nettoyer le système lancé automatiquement échoue	Exécutez le script : B03 – Nettoyer le système .

Autres problèmes

Problème	Script
État Erreur : une ou plusieurs barres de LED sont rouges et l'Vanquish User Interface indique que le système est dans l'état Erreur.	 Exécutez le script : B04 – Test automatique du système. Vérifiez si l'état général du système est OK (inême verte)
	 3. Si ce n'est pas le cas, vérifiez si le volet notification affiche des messages et résolvez le problème (reportez-vous à la section Messages () page 356)).
Les réglages du système ne correspondent pas	Pour régler le système ou des modules individuels du système aux valeurs par défaut d'usine, exécutez le script : E02 – Réinitialiser les réglages d'usine par défaut du système .
Il est impossible de récupérer les réglages du système et il n'a pas été correctement initialisé	Exécutez le script avec les paramètres appropriés à la configuration et au flux de travail : E01 – Initialiser le réglage du système.
La procédure de changement du circuit fluidique ou le flux de travail (A06 – Modifier le circuit fluidique / flux de tâches) n'ont pas été réalisés avec succès et seuls quelques scripts sont disponibles	

Recherchez les tableaux des sections suivantes qui correspondent au type de symptôme que vous observez, trouvez la cause possible et appliquez la solution proposée pour résoudre rapidement le problème.

8.3.1 Pics

Aucun pic

Cause possible	Mesure corrective
Appareil de préchauffage utilisé dans les applications LC nano ou capillaire	Retirez l'appareil de préchauffage (reportez-vous à la section Remplacement de l'appareil de préchauffage passif (▶ page 313)).
Mauvais capillaires utilisés entre la colonne et le détecteur ou entre le passeur d'échantillon et la colonne	Utilisez les capillaires adaptés à l'application (pour changer le flux de travail et les raccordements fluidiques, exécutez le script : A06 – Modifier le circuit fluidique / flux de tâches ou reportez-vous au <i>Schéma du débit</i>).
Présence de bulles d'air dans la boucle d'échantillonnage, l'aiguille ou le dispositif de mesure	Purgez le passeur d'échantillon.
Échantillon vide	Vérifiez le niveau d'échantillon dans l'échantillon. Si le niveau est trop bas, remplacez l'échantillon par un échantillon rempli.
Le solvant présent dans le réservoir de solvant ne correspond pas aux réglages concernant les solvants	Changez le solvant dans le réservoir de solvant ou changez les réglages concernant les solvants.

Dérive des pics

Cause possible	Mesure corrective
Température instable dans le compartiment à colonnes	 Vérifiez la température du compartiment à colonnes. Patientez jusqu'à ce que la température est stabilisée.
Après avoir changé de type de solvant, le solvant précédent n'a pas été complètement éliminé du passeur d'échantillon ou du système	 Si le script (B01 – Modifier les liquides / solvants) n'a pas complètement éliminé le solvant précédent, vérifiez si le bon paramètre a été configuré: Rafraîchir Uniquement = désactivé. Exécutez le script de nouveau : B01 – Modifier les liquides / solvants avec le paramètre : Rafraîchir Uniquement = désactivé.

8.3.1.1 Forme de pic

Traînée des pics

Cause possible	Mesure corrective
Volume hors colonne trop élevé	 Utilisez des raccords capillaires avec un diamètre interne adapté. Utilisez les capillaires Viper adaptés à l'application (reportez-vous au script : A06 – Modifier le circuit fluidique / flux de tâches).
Raccordements incorrects des capillaires	 Vérifiez le positionnement correct des raccords. Desserrez et resserrez les connecteurs Viper (reportez-vous à la section Connexion de capillaires nanoViper (▶ page 122)). Utilisez les capillaires Viper adaptés à l'application (reportez-vous au script : A06 – Modifier le circuit fluidique / flux de tâches). Remplacez le capillaire.
Volumes morts dans les raccordements des capillaires	 Exécutez le script : D02 – Tester le serrage du système (ce qui devrait être testé = l'échantillonneur). Inspectez le capillaire et remplacez-le, le cas échéant.

Élargissement de la bande

Cause possible	Mesure corrective
Volume hors colonne trop élevé	• Si les premiers pics sont plus larges que ceux qui éluent ultérieurement, vérifiez notamment le diamètre intérieur et la longueur du capillaire.
	 Utilisez un capillaire de plus faible volume pour raccorder le compartiment à colonnes au détecteur en aval.
Gradient de température dans la colonne	 Si vous utilisez des débits de > 50 μl/min, surtout avec des colonnes ID de 1 mm avec une gamme de débit entre 50 et 100 μl/min, utilisez un appareil de préchauffage.
	 SI un appareil de préchauffage est déjà installé, vérifiez qu'il l'est correctement. Au besoin, réinstallez-le (reportez-vous à la section Remplacement de l'appareil de préchauffage passif (> page 313)).
Capillaires	1. Exécutez le script : D01 – Tester la contre-pression du système.
partiellement obstrues	2. Remplacez les capillaires.
Boucle d'échantillonnage partiellement obstruée	Exécutez le script : D01 – Tester la contre-pression du système .
Volumes morts dans les raccordements des capillaires	1. Exécutez le script : D02 – Tester le serrage du système (ce qui devrait être testé = l'échantillonneur).
	2. Installez correctement les raccords des capillaires.
	3. Serrez les raccords des capillaires.
	4. Remplacez le capillaire.

Élargissement de la bande, temps mort accru

Cause possible	Mesure corrective
Volume après-colonne élevé	 Utilisez un capillaire avec les paramètres adaptés suivants : Diamètre intérieur Longueur Inspectez le capillaire. S'il est défectueux, remplacez-le.
Filtre du siège de l'aiguille obstrué	 1. Exécutez le script : D01 – Tester la contre-pression du système. 2. Remplacez le siège de l'aiguille, le cas échéant.

Pics triangulaires

Cause possible	Mesure corrective
Colonne surchargée (concentration d'échantillon trop élevée)	Diluez l'échantillon.

8.3.1.2 Pics fantômes, pics négatifs et pics parasites

Pics fantômes

Cause possible	Mesure corrective
Contamination (généralement : injecteur ou colonne)	Rincez le système, y compris la colonne, avec un solvant approprié (reportez-vous à la section Rinçage de la pompe en cas de contamination persistante uniquement () page 190).
Éluants dégradés, sales ou de mauvaise pureté/qualité	 Utilisez des solvants adaptés fraîchement préparés, par exemple, de qualité MS. Remplacez les réservoirs de solvant en utilisant des solvants adaptés fraîchement préparés (qualité MS) (exécutez le script : B01 – Modifier les liquides / solvants).

Pics négatifs

Cause possible	Mesure corrective
Composition différente du solvant d'échantillon et de la phase mobile	Dissolvez l'échantillon dans la phase mobile.
Air injecté/flacon vide	Vérifiez si le flacon est vide. Remplacez-le par un flacon rempli.

Pics parasites

Cause possible	Mesure corrective
Interférences électriques provenant d'autres modules du système	 Évitez l'emploi d'équipements générant de forts champs magnétiques ou électriques à proximité du système. Ne branchez pas d'appareils pouvant perturber l'alimentation électrique sur le même réseau que le système.
Température de colonne nettement supérieure au point d'ébullition de la phase mobile	En élaborant des méthodes, évitez que la température de colonne dépasse de manière significative le point d'ébullition de la phase mobile.
Solvant insuffisamment dégazé	Dégazez les solvants dans un bain à ultrasons.

8.3.1.3 Précision de l'aire des pics

Mauvaise précision de l'aire des pics/aires des pics trop petits

Cause possible	Mesure corrective
Présence de bulles d'air dans la boucle d'échantillonnage, l'aiguille ou le dispositif de mesure	Purgez le passeur d'échantillon.
Flacon vide	Vérifiez le niveau d'échantillon dans le flacon. Si le niveau est trop bas, remplacez le flacon par un flacon rempli.

Contamination ou transfert interéchantillons

Cause possible	Mesure corrective
Contamination présente dans le système	Rincez le système avec un solvant adapté. Si vous utilisez de l'eau comme solvant dans la voie A de la pompe, rincez-la avec de l'acétonitrile 80 %.
Transfert interéchantillons dans le passeur d'échantillon causé par la contamination de l'aiguille	 Vérifiez si l'une des conduites de rinçage de l'aiguille (étiquetées SWP et WWP) est pincée.
	 Vérifiez si l'un des réservoirs de rinçage de l'aiguille est vide. Si c'est le cas, remplacez-le en exécutant le script : B01 – Modifier les liquides / solvants. Utilisez un liquide de rinçage de l'aiguille adapté (reportez-vous à la section Remplissage des réservoirs de rinçage de l'aiguille (* page 131)).
	 Vérifiez si le rinçage de l'aiguille est désactivé dans le système de gestion de données chromatographiques. SI c'est le cas, activez le rinçage de l'aiguille.
	 Pour rincer l'aiguille et le siège de l'aiguille, exécutez le script : B03 – Nettoyer le système.
	5. Si le transfert interéchantillons se reproduit, remplacez l'unité d'aiguille et le siège de l'aiguille (reportez-vous à la section Remplacement du siège de l'aiguille et de l'unité de l'aiguille (l> page 254)).

Cause possible	Mesure corrective
Transfert interéchantillons dans le passeur d'échantillon causé par un siège de l'aiguille usé ou endommagé	Remplacez le siège de l'aiguille (reportez-vous à la section Remplacement du siège de l'aiguille et de l'unité de l'aiguille (▶ page 254)).

Analyse instable ou faible reproductibilité

Cause possible	Mesure corrective
Environnement instable	 Assurez-vous que la température et l'humidité de l'air sont constantes. Pour de l'information sur les conditions environnementales spécifiées, consultez le manuel d'utilisation du détecteur.
	Utilisez la thermostatisation des colonnes.
	 Si l'échantillon permet le refroidissement, refroidissez-le dans le passeur d'échantillon. Évitez les courants d'air.
Gradient non reproductible	 Vérifiez la valeur de %B. Si elle est à l'extérieur de la gamme 1-99, corrigez-la. Le système n'est pas correctement équilibré. Laissez-le s'équilibrer pendant 5 heures. Purgez la pompe.

Fuite

Cause possible	Mesure corrective
Raccords de capillaires incorrectement	 Exécutez le script : D02 – Tester le serrage du système (ce qui devrait être testé = le système).
installés ou insuffisamment serrés	Inspectez le capillaire et remplacez-le, le cas échéant.

Problèmes d'échantillons

Cause possible	Mesure corrective
Échantillon instable et se décompose	 Refroidissez l'échantillon dans le passeur d'échantillon. Utilisez un nouvel échantillon.
	 Changez les conditions environnementales.

Variation du volume d'injection

Cause possible	Mesure corrective
Aspiration de l'air du flacon par le passeur d'échantillon	La quantité d'échantillon dans le flacon est insuffisante ou la hauteur de l'aiguille est incorrecte.
	• Vérifiez le niveau d'échantillon dans le flacon. Si le niveau est trop bas, remplacez le flacon par un flacon rempli.
	 Pour vérifier si l'aiguille est obstruée, exécutez le script : D01 – Tester la contre- pression du système. Suivez les instructions.

Cause possible	Mesure corrective
Air emprisonné dans le trajet d'écoulement du passeur d'échantillon	Purgez le passeur d'échantillon.
Teneur en gaz de l'échantillon trop élevée ou saturée	Dégazez l'échantillon, si possible.
Aiguille de l'injecteur obstruée ou déformation de la pointe de l'aiguille	Exécutez le script : D01 – Tester la contre-pression du système . Suivez les instructions.
Fuite des joints de piston dans le dispositif de mesure du passeur d'échantillon	Exécutez le script : D02 – Tester le serrage du système (ce qui devrait être testé = l'échantillonneur). Suivez les instructions.
Fuite de la vanne d'injection ou d'autres composants du passeur d'échantillon	Exécutez le script : D02 – Tester le serrage du système (ce qui devrait être testé = l'échantillonneur). Suivez les instructions.
Siège de l'aiguille usé	Exécutez le script : D02 – Tester le serrage du système (ce qui devrait être testé = l'échantillonneur). Suivez les instructions.

8.3.2 Pression

Pulsation de pression ou pression non constante

Cause possible	Mesure corrective
Système encrassé par des particules	Les échantillons et une pureté insuffisante des solvants sont souvent la cause de l'encrassement du système par des particules. Dans les applications LC nano et LC capillaire, les capillaires de diamètres encore plus petits nécessitent des solvants et des échantillons encore plus purs.
	 Applications LC nano et LC capillaire : Pour éviter la contamination de l'échantillon, procédez comme suit :
	 Utilisez les échantillons avec du septum sans particules du kit d'expédition ou ceux énumérés dans la Recommandations sur les bouchons de flacons et les couvercles des plaques à puits (> page 167).
	 Réduisez le nombre de répétitions d'un échantillon bouché à < 10.
	 Rincez l'aiguille et/ou intensifiez le rinçage de l'aiguille dans le système de gestion de données chromatographiques.
	 Exécutez le script : D01 – Tester la contre-pression du système. Suivez les instructions.
Solvant dégradé	Utilisez du solvant fraîchement préparé.

Contre-pression élevée

Cause possible	Mesure corrective
Encrassement ou endommagement par torsion d'un ou plusieurs capillaires du système	Exécutez le script : D01 – Tester la contre-pression du système . Suivez les instructions.
Colonne contaminée ou encrassée	 Exécutez le script : D01 – Tester la contre-pression du système. Suivez les instructions.
	2. Si le test réussi, la colonne ou le capillaire de l'appareil en aval est encrassé.
	3. Remplacez la colonne ou le capillaire vers l'appareil en aval.
Encrassement de la	1. Inspectez la vanne à la recherche de signes de colmatage.
vanne de commutation pour colonne	 Remplacez la vanne, le cas échéant (reportez-vous à la section Remplacement de la vanne de commutation Vanquish (> page 305) ou Remplacement de la vanne de commutation VICI (> page 308)).
Encrassement de l'appareil de préchauffage	Remplacez l'appareil de préchauffage (reportez-vous à la section Remplacement de l'appareil de préchauffage passif () page 313)).
Contamination présente dans le système	Rincez le système en purgeant la pompe et le passeur d'échantillon avec un solvant adapté.
Le siège de l'aiguille obstrué	Si le script de test de la contre-pression affiche un message indiquant qu'une grande contre-pression est détectée au niveau du siège de l'aiguille, exécutez le script : C21 – Nettoyer ou remplacer le siège et l'unité d'aiguille avec le paramètre : Nettoyer le siège de l'aiguille .
Le solvant présent dans le réservoir de solvant ne correspond pas aux réglages concernant les solvants	Changez le solvant dans le réservoir de solvant ou changez les réglages concernant les solvants.

Contre-pression basse

Cause possible	Mesure corrective
Fuite dans le système	Exécutez le script : D02 – Tester le serrage du système (ce qui devrait être testé = le système). Suivez les instructions.
Air emprisonné dans le passeur d'échantillon	Exécutez le script : CO4 – Purger l'échantillonneur.
Le solvant présent dans le réservoir de solvant ne correspond pas aux réglages concernant les solvants	Changez le solvant dans le réservoir de solvant ou changez les réglages concernant les solvants.

8.3.3 Variation des temps de rétention

Réduction des temps de rétention

Cause possible	Mesure corrective
Composition de solvant inappropriée	 Vérifiez l'âge du solvant prémélangé. Purgez la pompe. En préparant l'utilisation du système, laissez-le s'équilibrer pendant 5 à 24 heures. Exécutez le script : C01 – Régler les décalages du capteur de débit de la pompe.

Augmentation des temps de rétention

Cause possible	Mesure corrective
Composition de solvant inappropriée	 Vérifiez l'âge du solvant prémélangé. Purgez la pompe. En préparant l'utilisation du système, laissez-le s'équilibrer pendant 5 à 24 heures. Exécutez le script : C01 – Régler les décalages du capteur de débit de la pompe.
Débit décroissant	 Exécutez le script : D02 – Tester le serrage du système (ce qui devrait être testé = le système). Suivez les instructions. Vérifiez les paramètres de débit.

Dispersion des temps de rétention

Cause possible	Mesure corrective
Fluctuations de pression de la pompe	Exécutez le script : D02 – Tester le serrage du système (ce qui devrait être testé = le système).
Air emprisonné dans le passeur d'échantillon	Exécutez le script : CO4 – Purger l'échantillonneur avec le paramètre : Liquide de rinçage de l'aiguille faible pour rinçage interne.

Décalages des temps de rétention pendant une période prolongée

Cause possible	Mesure corrective
Après avoir changé les solvants, le solvant précédent n'a pas été complètement éliminé de la pompe.	Purgez la pompe.
Température de colonne, d'échantillon ou de solvant instable	Patientez jusqu'à ce que la température de la colonne, de l'échantillon ou du solvant est complètement équilibrée.

Cause possible	Mesure corrective
Il est possible que les capteurs de débit aient une petite dérive de la ligne de base ou décalage.	Exécutez le script : CO1 – Régler les décalages du capteur de débit de la pompe.
Les solvants prémélangés sont trop vieux.	 Utilisez un éluant fraîchement préparé. Utilisez des solvants adaptés fraîchement préparés (par exemple, de qualité MS). Purgez la pompe.
Appareil de préchauffage utilisé dans les applications LC nano	Retirez l'appareil de préchauffage (reportez-vous à la section Remplacement de l'appareil de préchauffage passif () page 313)).
Volume après-colonne erroné	Vérifiez le volume du capillaire connecté à l'orifice de SORTIE de la colonne.
Le variant de la boucle d'échantillonnage ne correspond pas au varient configuré	Vérifiez si le réglage de la boucle d'échantillonnage correspond au réglage de la boucle installée et corrigez-le, le cas échéant.
Le solvant présent dans le réservoir de solvant ne correspond pas aux réglages concernant les solvants	Changez le solvant dans le réservoir de solvant ou changez les réglages concernant les solvants.

Temps de rétention à l'extérieur de la gamme attendue

8.3.4 Ligne de base

Dérive importante de la ligne de base

Cause possible	Mesure corrective
Système insuffisamment équilibré	 Rincez le système jusqu'à équilibrage. Habituellement, un volume de 5 à 10 fois celui de la colonne est suffisant.
	 Si le rinçage du système ne résout pas le problème, purgez la pompe et laissez le système s'équilibrer pendant au moins 5 heures.
Environnement instable	Assurez-vous que la température et l'humidité de l'air sont constantes.
L'absorption de l'éluant varie au cours du gradient	Les additifs absorbants peuvent modifier le spectre d'absorption en fonction du solvant. Envisagez l'utilisation de concentrations en additifs variables pour limiter la dérive.
Contamination du passeur d'échantillon	Exécutez le script : B03 – Nettoyer le système .
Contamination de la colonne	Nettoyez ou remplacez la colonne.

Cause possible	Mesure corrective
Vieux solvant présent dans la tête du dispositif de mesure ou la boucle d'échantillonnage	Pour purger le dispositif de mesure ou la boucle d'échantillonnage, exécutez le script : C04 – Purger l'échantillonneur avec le paramètre Liquide de rinçage de l'aiguille fort pour rinçage interne .
	En général, lorsque vous remplacez les réservoirs de solvant, exécutez le script : B01 – Modifier les liquides / solvants.
Solvants dégradés ou non homogènes	 Avant de commencer une analyse, assurez-vous que les solvants sont déjà homogénéisés dans les réservoirs.
	 Dans les solvants aqueux, de microorganismes peuvent se développer et dégrader le solvant. Pour remplacer le solvant par un nouveau réservoir de solvant contenant du solvant fraîchement préparé, exécutez le script : B01 – Modifier les liquides / solvants. Suivez les instructions.

Fluctuations non périodiques de la ligne de base, niveau de bruit élevé

Cause possible	Mesure corrective
Air emprisonné dans le système	Purgez la pompe et le passeur d'échantillon.
Solvant dégradé ou de mauvaise qualité	 Utilisez un éluant fraîchement préparé. Utilisez des solvants adaptés (de qualité MS). Pour changer de solvant, exécutez les scripts : A01 – Définir le type de solvant de la pompe, puis B01 – Modifier les liquides / solvants.

Fluctuations périodiques de la ligne de base, pulsation

Cause possible	Mesure corrective
Fluctuations de pression de la pompe	 Purgez la pompe. Exécutez le script : D01 – Tester la contre-pression du système.
Air emprisonné dans le système	Purgez la pompe.

8.3.5 Temps de chargement de la colonne de piégeage prolongé

Cause possible	Mesure corrective
Colonne de piégeage vieillie ou obstruée	1. Exécutez le script : D01 – Tester la contre-pression du système avec le paramètre Tester uniquement la résistance de la colonne de piégeage.
	2. Écrivez la valeur de résistance et comparez-la à un ancien résultat.
	3. Si la résistance est trop élevée pour votre application, effectuez le test de contre- pression complet. Exécutez le script : D01 – Tester la contre-pression du système .
	4. Vérifiez le résultat du test et remplacez le composant obstrué, le cas échéant.
	 Si vous avez remplacé un composant, effectuez à nouveau le test de contre- pression : D01 – Tester la contre-pression du système.
Le siège de l'aiguille obstrué	Reportez-vous à la sous-section <i>Contre-pression élevée</i> de la section Pression () page 386).

8.3.6 Régulation de la température

8.3.6.1 Problèmes de la régulation de la température du compartiment à échantillons

La température du compartiment à colonnes ne change pas alors que la valeur de consigne n'est pas atteinte

Cause possible	Mesure corrective	
Température ambiante trop élevée ou différence entre la température de consigne et la température ambiante inférieure à celle précisée dans les caractéristiques techniques	 Abaissez la température ambiante (en climatisant le local par exemple). Sélectionnez une valeur de consigne plus élevée pour la température. 	
Régulation de la température désactivée	Activez la régulation de la température.	
Obstruction de l'ouverture de ventilation	 Ouvrez les portes du compartiment à colonnes et patientez jusqu'à l'arrêt des ventilateurs. Vérifiez si les ouvertures de ventilation du compartiment à colonnes sont obstruées. Retirez toute particule obstruante des ouvertures de ventilation. 	
Entrée de l'air ambiant dans la chambre à colonnes	 Une mauvaise étanchéité de la chambre à colonnes peut nuire au chauffage et au refroidissement du compartiment à colonnes et provoquer une très forte condensation. 1. Vérifiez si les portes de devant ferment correctement. 2. Vérifiez si les portes sont bien étanches. 3. Vérifiez si le joint de la porte est endommagé. Si c'est le cas, remplacez la porte (reportez-vous à la section Remplacement des portes () page 333)). 	
Compartiment à colonnes défectueux	Prenez contact avec l'assistance technique.	

8.3.6.2 Problèmes de la régulation de la température du passeur d'échantillon

Le	passeur	d'échantillon	n'atteint pas	la température	de consiane
LC	pusseur	a centantinon	n attent pas	ia temperature	ac consigne

Cause possible	Mesure corrective
Température ambiante ou humidité trop élevée	 Vérifiez si le couvercle isolant est installé. Abaissez la température ambiante (en ventilant le local par exemple) et/ou l'humidité ambiante.

8.3.7 Alimentation et communications (modules du système)

Les problèmes et les actions correctives décrits dans cette section portent sur les modules du système (par exemple, le socle du système, la pompe, le passeur d'échantillon et le compartiment à colonnes [le cas échéant]). Pour obtenir des informations sur les problèmes d'alimentation et de communications du contrôleur du système et de l'interface utilisateur, reportez-vous à la section Problèmes en lien avec le contrôleur du système, le dispositif d'affichage et l'interface utilisateur () page 393).

Aucune partie du système n'est alimentée (toutes les barres de LED sont éteintes [sombres])

Cause possible	Mesure corrective
Bouton d'alimentation du système en position Arrêt	Mettez le système sous tension au moyen du bouton d'alimentation situé sur le devant gauche du socle du système.

La LED d'état du contrôleur du système est éteinte (sombre)

Cause possible	Mesure corrective	
Bouton d'alimentation du contrôleur du système désactivé	Mettez le contrôleur du système sous tension au moyen du bouton d'alimentation supérieur situé sur le socle du système.	
Interrupteur d'alimentation principal du contrôleur du système désactivé	Mettez le contrôleur du système sous tension au moyen de son interrupteur d'alimentation principal.	
Module non branché au secteur	Branchez le cordon d'alimentation.	
Erreur du système électronique	Prenez contact avec l'assistance technique.	

Aucune alimentation en électricité du module (barre de LED d'état éteinte [sombre])

Cause possible	Mesure corrective	
Alimentation du module éteinte	Mettez le module sous tension au moyen de son interrupteur d'alimentation principal.	
Module non branché au secteur	Branchez le cordon d'alimentation.	
Fusible grillé	Remplacez le fusible (reportez-vous à la section Remplacement des fusibles de l'alimentation principale () page 332)).	
Erreur du système électronique	Prenez contact avec l'assistance technique.	

Impossible de mettre hors tension un ou plusieurs modules avec le bouton d'alimentation sur le socle du système

Cause possible	Mesure corrective
Câbles System Interlink manquants ou défectueux	Vérifiez les branchements de System Interlink entre le socle du système et les modules, reportez-vous à la section Connexion de câbles System Interlink () page 113).

Impossible de piloter le module à partir du système de gestion de données chromatographiques

ions USB entre le système et l'ordinateur, reportez-vous à la de câbles USB (▶ page 108). Si le module utilise un situé sur un autre module, assurez-vous que le module avec sous tension. on USB sur l'ordinateur. Une connexion USB standard 2.0 est

8.3.8 Problèmes en lien avec le contrôleur du système, le dispositif d'affichage et l'interface utilisateur

Symptôme	Cause possible	Mesures correctives
Clignotement du dispositif d'affichage	Dispositif d'affichage mal connecté	Vérifiez si le connecteur du câble du dispositif d'affichage est serré manuellement au port Display Data du contrôleur du système.
Dispositif d'affichage non fonctionnel	Dispositif d'affichage non connecté au contrôleur du système	 Vérifiez si le dispositif d'affichage est connecté au contrôleur du système. Si ce n'est pas le cas, passez à l'étape suivante. Mettez le contrôleur du système hors tension
		 3. Connectez le dispositif d'affichage au contrôleur du système (reportez-vous à la section Connexion des câbles du dispositif d'affichage (> page 111)). 4. Mettez le contrôleur du système sous tension.
Le contrôleur du système ne peut pas être désactivé avec son bouton d'alimentation	Câble VSC court non branché	Assurez-vous que le câble VSC court relie le port VSC du socle du système au port VSC du contrôleur du système.
Le contrôleur du système est sous tension, mais la LED d'état est éteinte (sombre)		

Symptôme	Cause possible	Mesures correctives
Le contrôleur du système ne peut pas être activé alors que vous avez appuyé sur le bouton d'alimentation principal		
Aucun script disponible	-	Mettez le contrôleur du système hors tension, patientez pendant 5 secondes et allumez-le à nouveau.
Aucun module du système n'apparaît sur l'Vanquish User Interface ou il y a des scripts manquant dans l'Vanquish User Interface	Système empilé non connecté au port System Interlink du contrôleur du système	Connectez le système empilé au port System Interlink sur le panneau court du contrôleur du système (voir Connexion de câbles System Interlink () page 113)).
	Système empilé non connecté au bon port System Interlink du contrôleur du système	Connectez le système empilé au port System Interlink sur le panneau court du contrôleur du système (voir Connexion de câbles System Interlink () page 113)).
Le module du système unique n'apparaît pas sur l'Vanquish User Interface ou il y a des scripts manquant dans l'Vanquish User Interface	Module du système unique non connecté à l'interface System Interlink	 Vérifiez si le module système manquant est connecté à l'interface System Interlink du système. Si ce n'est pas le cas, intégrez-le dans le bus System Interlink du système empilé (reportez-vous à la section Connexion de câbles System Interlink () page 113)). Vérifiez si le module du système est mis sous tension au moyen de son interrupteur d'alimentation principal. Remplacez le câble System Interlink.
	Version du microprogramme du module et / ou du contrôleur du système incorrecte ou incohérence des versions de microprogramme	Installez les versions correctes et correspondantes du microprogramme sur le module et / ou sur le contrôleur du système. Prenez contact avec l'assistance technique de Thermo Fisher Scientific pour obtenir de l'aide.
Le numéro de série sur l'Vanquish User Interface est Indéfini	Le numéro de série n'a pas été défini après le remplacement d'un module à glissières par un nouveau.	Faites programmer le numéro de série du système dans le module du système par un technicien d'entretien Thermo Fisher Scientific.

8.4 Recherche de fuites de liquide dans le système

Vérifiez régulièrement si des fuites des liquides suivants sont présentes dans le système :

Liquide	Suivez les instructions de
Solvant	Cette section
Liquide de rinçage des joints	Test d'étanchéité du système de rinçage des joints () page 282)
Solvant et liquide de rinçage de l'aiguille des vannes du passeur d'échantillon	Inspection des vannes du passeur d'échantillon à la recherche de fuites (▶ page 301)
Eau de condensation provenant du passeur d'échantillon	Test des tubes de la pompe d'évacuation (▶ page 268)

À quel moment exécuter cette opération ?

- Régulièrement (comme partie du plan de maintenance) :
 - Après le remplacement des réservoirs de solvant
 - Après le changement du flux de travail
- Un capteur de fuites signale une fuite.
- Problèmes de chromatographie :
 - Pics trop petits
 - Variation dans les temps de rétention

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 179).

Procédez comme suit

1. Dépend des éléments suivants :

À quel moment exécuter cette opération ?	Action
Le test est régulièrement effectué (dans le cadre du plan de maintenance)	Exécutez le script : D02 – Tester le serrage du système (ce qui devrait être testé = le système) .
Le test est effectué, car : • un capteur de fuites signale une fuite ; • des problèmes de chromatographie sont observés.	 Exécutez le script : D02 – Tester le serrage du système (ce qui devrait être testé = le système). Exécutez le script : E04 – Test de fuite détaillé (pompe)

- 2. Contrôlez le résultat du test. Si le script signale une fuite, réparez-la en suivant les instructions à l'écran.
- 3. Pour vérifier que la fuite est complètement réparée, exécutez le script de nouveau et vérifiez les résultats.
9 Caractéristiques techniques

Le présent 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 du système.

9.1 Caractéristiques physiques

9.1.1 Spécifications environnementales

Les conditions environnementales sont spécifiées comme suit :

Туре	Caractéristique technique
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
Catégorie de surtension	Ш

9.1.2 Exigences relatives à l'alimentation électrique

Les exigences relatives à l'alimentation électrique sont spécifiées comme suit :

Produit	Exigences relatives à l'alimentation électrique
Pompe	100 – 240 V CA, ± 10 % ; 50/60 Hz ; max. 525 W / 550 VA
Passeur d'échantillon	100 – 240 V c.a., ± 10 % ; 50/60 Hz ; max. 525 W / 550 VA
Compartiment à colonnes	100 – 240 V c.a., ± 10 % ; 50/60 Hz ; max. 525 W / 550 VA
Contrôleur du système	100 – 240 V CA, ± 10 % ; 50 /60 Hz, max. 50 W / 150 VA

9.1.3 Dimensions

Les dimensions du système et des modules du système sont spécifiées comme suit :

Produit	Dimension (hauteur × largeur × profondeur)	
Système	Configuration de base ¹ : 74 x 45 x 65 cm (dispositif d'affichage installé sur le passeur d'échantillon et placé devant le bac à solvants)	Configuration de base et compartiment à colonnes : 82 x 45 x 65 cm (dispositif d'affichage installé sur le passeur d'échantillon et placé devant le compartiment à colonnes)
Pompe	19,2 x 42 x 62 cm	
Passeur d'échantillon	29 x 42 x 62 cm	
Compartiment à colonnes	Dimensions extérieures	15,9 x 42 x 62 cm
	Chambre à colonnes	7,8 x 34,3 x 3,7 cm

¹ : La configuration de base comprend le socle du système avec le contrôleur du système, la pompe, le passeur d'échantillon et le dispositif d'affichage

9.1.4 Niveau de pression acoustique

Les niveaux de pression acoustique sont spécifiés comme suit :

Produit	Niveau de pression acoustique
Système	< 75 dB(A), généralement < 55 dB(A)
Pompe	< 70 dB(A), généralement < 50 dB(A)
Passeur d'échantillon	< 70 dB(A), généralement < 50 dB(A)
Compartiment à colonnes	< 42 dB(A), généralement < 40 dB(A)

9.1.5 Poids

Les poids du système et des modules du système dont spécifiés comme suit :

Produit	Poids	
Système	Configuration de base ¹ : Env. 66 kg	Configuration de base et compartiment à colonnes : Env. 79 kg
Pompe	24 kg	
Passeur d'échantillon	26 kg	
Compartiment à colonnes	Sans vannes : 10,7 kg ; avec 2 vannes Vanquish : 11,6 kg ; avec 2 vannes VICI : 13,1 kg	
¹ : La configuration de base comprend le socle du système avec le contrôleur du système, la pompe, le passeur d'échantillon et le dispositif d'affichage		

9.2 Caractéristiques de performance

9.2.1 Caractéristiques de performance du système

La performance du système est spécifiée comme suit :

Туре	Caractéristique technique
Communication	Le socle du système est muni des ports suivants : 1 port System Interlink (connecteur RJ45-8) 1 port VSC (connecteur RJ-12) Pour les ports des modules du système, reportez-vous aux sections sur les caractéristiques de performance des modules du système.
Commande	Vanquish User Interface Panneau de commande doté de boutons permettant d'exécuter certaines fonctions depuis le module du système. Le système peut être piloté avec un système de gestion de données chromatographiques. Pour en savoir plus, veuillez prendre contact avec le service commercial de Thermo Fisher Scientific.
Matériaux du trajet d'écoulement	Pour les matériaux du trajet d'écoulement, reportez-vous aux sections sur les caractéristiques de performance de la pompe, du passeur d'échantillon et du compartiment à colonnes.
Informations sur les solvants et les additifs	Reportez-vous à la section Utilisation de solvants et d'additifs () page 145).
Biocompatibilité	Oui
Fonctions de sécurité	Détection de fuites et gestion sécurisée des fuites, surveillance de pression excessive
Fonctions de bonnes pratiques de laboratoire (BPL)	Les fonctions Predictive Performance permettent de programmer les opérations de maintenance en fonction des conditions réelles de fonctionnement et d'utilisation des modules du système. Tous les paramètres du système sont consignés dans le volet notification.

9.2.2 Caractéristiques de performance de la pompe

Туре	Spécifications
Principe de fonctionnement	Pompe à double piston en série avec entraînements de piston indépendants
Compensation de la compressibilité	Automatisation complète, indépendante de la composition de la phase mobile
Plage de débit (réglable)	1 nL/min – 100 μL/min par incréments de 1 nL/min
Exactitude du débit	Généralement : ± 3 %
	Condition : 0,1 à 100 μL/min pour l'un des 4 solvants pré-étalonnés, mesuré à une pression de système de 20 MPa
Déviation entre les modules	< 5 % pour les débits de > 300 nL/min – 10 µL/min
	Conditions : mesurée en différences de temps de rétention de peptides de protéines digestes TGQAPGFSYTDANK cytochrome C et un gradient de 30 min.
Précision du temps de rétention	\leq 0,2 % RSD ou \leq 0,1 SD min, le plus élevé des deux
	Conditions : pour les peptides de protéines digestes TGQAPGFSYTDANK cytochrome C pour un gradient de 30 min à 90 min.
Plage de pression	Pression de la pompe : 2 – 150 MPa (20 – 1500 bar, 300 – 21750 psi)
	Pression de pompe de 150 MPa jusqu'à 1 μL/min. Avec un débit supérieur à 1 μL/min, la pression maximale dépend du solvant et diminue de manière linéaire jusqu'à 70 MPa à 100 μL/min pour de l'acétonitrile 80 % sur la voie B.
Nombre de voies de solvants	2
Volume de course maximal	96 μL
Volume de délai (contribution de la pompe au volume de retard des gradients du système)	25 nL
Dégazage	2 voies pour les liquides de rinçage de l'aiguille du passeur d'échantillon
Communication	USB :
	1 port USB (USB 2.0, connecteur de type « B »)
	1 concentrateur USB muni de 3 ports (USB 2.0, connecteurs de type « A »)
	Interface E/S :
	2 ports Dig I/O (mini DIN), chacun fournissant une entrée et une sortie relais
	System Interlink :
	2 ports System Interlink (connecteurs RJ45-8)
Matériaux dans le trajet d'écoulement analytique	Céramique, DLC (carbone de type diamant), fluoropolymères, MP35N, PEEK, titane, PE-UHMW
	Pour de l'information sur la résistance chimique des matériaux, consultez la littérature technique.

9.2.3 Caractéristiques de performance du passeur d'échantillon

Туре	Caractéristique techni	que
Méthode d'injection	Avec division du flux, injection directe ou emprisonner et éluer	
Nombre d'unités d'injection	1	
Plage de pression	5 – 151 MPa (50 – 1 51	17 bars, 700 – 22 002 psi)
Plage de volume d'injection	Boucle d'échantillonnage de 25 μl (standard)	0,01 à 25 $\mu l,$ incréments de 0,01 μl minimum
	Boucle d'échantillonnage de 100 μl (étendue)	0,01 à 100 μl, incréments de 0,01 μl minimum Grande quantité de piégeage grâce à multiples cycles de prélèvement et de cycles de piégeage : jusqu'à 500 μl
Exactitude du volume d'injection	Typiquement : ± 0,5 %	pour un volume d'injection de 5 μ l et de l'eau
Précision du volume d'injection	0,05 à 0,20 μl	<5,0 %
	0,20 à 0,50 μl	<1,5 %
	0,50 à 2,00 μl	<0,5 %
	Supérieure à 2 µl	<0,25 %
Caractéristiques de faible perte d'échantillon	2,5 μl sur 3 μl avec des flacons adaptés (reportez-vous à la section Recommandations sur les bouchons de flacons et les couvercles des plaques à puits () page 167))	
Linéarité	r > 0,9999 (caféine dans de l'eau)	
Rinçage de l'aiguille	Trempage et rinçage en continu de la surface externe de l'aiguille avec 2 liquides de rinçage de l'aiguille indépendants.	
	Rinçage de la boucle d'échantillonnage et de la surface intérieure de l'aiguille avec du liquide de rinçage de l'aiguille fort de l'orifice de rinçage (SWP). Ensuite, rinçage de la boucle d'échantillonnage et de la surface intérieure de l'aiguille avec du liquide de rinçage de l'aiguille faible du dispositif de mesure (W).	
Nombre de liquides de rinçage de l'aiguille	2	
Volume de délai	Pour un volume de boucle d'échantillonnage de 25 μl :	
(contribution du passeur d'échantillon au volume de retard des gradients du système)	En position d'injection	55 μl
and Bradients an systemer	En position bypass	51 nl
Transfert interéchantillons	< 0,002 % avec la caféi	ne
Lecteur de code-barres	Oui	
Lecture de codes-barres des fonctions d'automatisation	Lecture de codes-barres : • Détection des segments vides • Vérification des portoirs / plaques à puits • Gestion d'inventaire	

La performance du passeur d'échantillon est spécifiée comme suit :

Туре	Caractéristique technique
Thermostatisation du compartiment à échantillons : plage de température	De 4 à 40°C, refroidissement ≥ 23 K au-dessous de la température ambiante avec une humidité relative < 80 %
Exactitude de la température	-2°C et +4°C
Stabilité de la température	±1°C
Capacité d'accueil d'échantillons	4 segments disponibles pour les portoirs d'échantillons ou plaques à puits au format SBS ; plus des porte-flacons supplémentaires pouvant accueillir des flacons d'une capacité de (nombre de flacons x D.E.) : 12 x 22,5 mm
Capacité d'accueil de flacons	• 54 flacons d'un D.E. de 12 mm
	• 96 flacons d'un D.E. de 6 mm
	• 96 flacons d'un D.E. de 7 mm
	• 96 flacons d'un D.E. de 8 mm
	• 16 flacons d'un D.E. de 15 mm
Capacité d'accueil de plaques à puits	Plaques 96 et 384 puits au format SBS
Volume d'échantillon minimum requis	2 μl pour un volume d'injection de 1 μl
Communication USB	1 port USB (USB 2.0, connecteur de type "B")
	1 concentrateur USB équipé de 3 ports (USB 2.0, connecteurs de type "A")
Interface E/S	2 ports E/S numériques (mini DIN), chacun fournissant une entrée numérique et une sortie relais
System Interlink	2 ports System Interlink (connecteurs RJ45-8)
Matériaux dans le trajet	Titane, céramiques, PEEK, MP35N, DLC, saphir, fluoropolymères
d'écoulement analytique	Avis : pour obtenir des informations sur le liquide de rinçage des joints à utiliser, reportez-vous à la section Remplissage du réservoir de liquide de rinçage des joints (▶ page 134).
	Pour obtenir des informations sur la résistance chimique des matériaux, consultez la littérature technique.
Matériaux dans le trajet d'écoulement de rinçage de l'aiguille	PP (polypropylène), PE, FFPM (perfluoroélastomère), FFKM, PEEK, PA (polyamides), PK (polycétone), TPE (élastomère thermoplastique)
Matériaux dans le trajet d'écoulement de rinçage des joints	Titane, céramique, PEEK, UHMW, PE, PP, PTFE, silicone

9.2.4 Caractéristiques de performance du contrôleur du système

La performance du contrôleur du système est spécifiée comme suit :

Туре	Caractéristique technique
Interfaces E/S	1 port LAN
	2 ports USB (désactivés)
	4 ports System Interlink (2 sont désactivés)
System Interlink	2 ports System Interlink (1 paire, connecteurs RJ45-8)

9.2.5 Caractéristiques de performance du compartiment à colonnes

Туре	Spécifications
Plage de pression	Vanne Vanquish : 50 – 1 500 bar Vanne VICI C82 : 50 – 1 034 bar Vanne VICI C2, C2N : 50 – 345 bar
Plage de température	Min. : 5 K au-dessus de la température ambiante Max. : 80°C sans vannes installées Max. : 60°C avec des vannes Vanquish installées Max. : 50°C avec des vannes VICI installées
Exactitude de la température	± 0,5°C
Stabilité de la température	± 0,05°C
Précision de température	± 0,1°C
Délai de chauffage	De 35°C à 65°C ± 1°C en moins de 12 minutes
Vannes de commutation pour colonne	Jusqu'à deux vannes : Vanne Vanquish 2 positions 6 ports, 1 500 bar Vanne VICI C82 2 positions 6 ports, 1 034 bar Vanne VICI C2 2 positions 6 ports, 345 bar Vanne VICI C2N 2 positions 6 ports, 345 bar Pour des information supplémentaires sur les vannes de commutation pour colonne offertes, reportez-vous à la section Accessoires en option (▶ page 413).
Communication USB	1 port USB (USB 2.0, connecteur de type "B") 1 concentrateur USB muni de 3 ports (USB 2.0, connecteurs de type "A")
Interface E/S	2 ports Dig I/O (mini DIN), chacun fournissant une entrée numérique et une sortie relais
System Interlink	2 ports System Interlink (connecteurs RJ45-8)
Matériaux du trajet d'écoulement	Vanne Vanquish : céramique, DLC, titanium Vanne VICI C82 : acier inoxydable avec revêtement inerte, composite à base de PAEK Vanne VICI C2, C2N : PAEK, composite à base de PAEK Capillaires : MP35N (pour en savoir plus, consultez les spécifications de l'appareil de préchauffage) Pour obtenir des informations sur la résistance chimique des matériaux, consultez la documentation technique.

La performance du compartiment à colonnes est spécifiée comme suit :

Appareil de préchauffage passif

Туре	Spécifications
Plage de température	Identique à la chambre à colonnes
Matériaux dans le trajet d'écoulement	MP35N
Biocompatibilité	Oui
Volume chauffé	1 μL
Volume total	6 µL

10 Accessoires, consommables et pièces de rechange

Le présent chapitre répertorie les accessoires par défaut expédiés avec le système et les accessoires disponibles en option. Il fournit également des renseignements sur le réapprovisionnement en consommables et en pièces de rechange.

10.1 Généralités

Le système 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

Cette section liste le contenu du kit d'expédition. 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 du système.

Pour obtenir les informations de réapprovisionnement, reportez-vous à la section Consommables et pièces de rechange () page 417).

10.2.1 Système

Capillaires

Pour de l'information détaillée concernant l'utilisation des capillaires, reportez-vous au schéma de flux du kit d'expédition.

	Élément	Quantité dans l'expédition
-	Capillaire, nanoViper, gainage en verre silice fondue/PEEK, 1/32 po, D.I. x D.E. x longueur 20 μm x 360 μm x 550 mm, pour les applications nano LC	2
	Capillaire colonne – détecteur, nanoViper, gainage en silice fondue/PEEK, 1/32 po, D.I. x D.E. x longueur 20 μm x 280 μm x 1 m, pour les applications nano LC	1
	Capillaire colonne – détecteur, nanoViper, gainage en silice fondue/PEEK, 1/32 po, D.I. x D.E. x longueur 50 μm x 360 μm x 550 mm, pour les applications capillaires et micro LC	4
	Capillaire pompe – passeur d'échantillon, nanoViper, gainage en silice fondue/PEEK, 1/32 po, D.I. x D.E. x longueur 50 μm x 360 μm x 350 mm, pour les applications capillaires et micro LC	1
	Capillaire passeur d'échantillon – colonne/compartiment à colonnes, nanoViper, gainage en silice fondue/PEEK, 1/32 po, D.I. x D.E. x longueur 50 μm x 280 μm x 1 mm, pour les applications capillaires et micro LC	1
	Capillaire colonne – détecteur, silice fondue, D.I. x D.E. x longueur 20 μm x 280 μm x 1,5 m, pour les applications nano LC	1
	Capillaire colonne – détecteur, silice fondue, D.I. x D.E. x longueur 50 μm x 280 μm x 1,5 m, pour les applications capillaires et micro LC	1

Accessoires des capillaires

Élément	Quantité dans l'expédition
Raccord-union Viper	1
Bouchon de raccordement, Viper, titane	3
Molette pour raccords nanoViper	2

Accessoires du système de rinçage

Élément	Quantité dans l'expédition
Tubes, silicone	3 m
Raccord de tube, droit (paquet de 5)	1

Réservoirs de liquide et accessoires des réservoirs de liquide

Élément	Quantité dans l'expédition
Insertion de bouteille	1
Réservoir, 1 L, avec bouchon	5
Réservoir, 0,25 L, avec bouchon	4
Bouchon de réservoir	10
 Guide de maintien (diamètre de 1,6 mm) pour garder les conduites suivantes dans le bouchon de réservoir : Conduites de solvant (entre le réservoir et la pompe) Conduite de liquide de rinçage de l'aiguille (entre le réservoir et le dégazeur du liquide de rinçage de l'aiguille dans la pompe et entre le réservoir et la pompe de rinçage du passeur d'échantillon) 	10
Guide de maintien (diamètre de 3,0mm) pour garder la conduite de rinçage des joints dans le bouchon de réservoir	10
Capuchon protecteur pour refermer les trous ouverts dans le bouchon de réservoir (paquet de 20)	2
Filtre de conduite de solvant, porte-filtre (fritté ou adaptateurs de conduite de solvant non compris)	6
Adaptateur de conduite de solvant	6
Filtre de conduite de solvant, fritté de filtre, 10 μm	10

Outils

Élément	Quantité dans l'expédition
Outil de manipulation des joints	1
Seringue, plastique, 12 mL	2
Tournevis dynamométrique, Viper, pièce de base avec adaptateur d'extension et embouts pour serrer les capillaires nanoViper et le siège de l'aiguille du passeur d'échantillon	1
Coupe-tube pour les tubes PEEK	1
Coupe-tube pour les tubes en silice fondue	1

Câbles

Élément	Quantité dans l'expédition
Câble de signalement numérique d'entrée et de sortie, 5 m	1
Câble LAN (RJ45), 5 m	1
Câble USB, de type A à type B, USB 2.0 à haut débit, longueur 0,5 m	1
Pour connecter la pompe au passeur d'échantillon.	
Câble USB, de type A à type B, USB 2.0 à haut débit, longueur 5 m Pour connecter la pompe à l'ordinateur.	1

Éléments spécifiques au passeur d'échantillon

Élément	Quantité dans l'expédition
Siège de l'aiguille	1
Portoir d'échantillons, accueillant 54 échantillons d'un D.E. de 12 mm	4
Remarque : les portoirs d'échantillons sont munis d'un code- barres 2D pour l'identification du type de portoir Vanquish.	
Kit de récipients d'échantillonnage, avec plaques à puits, tapis, ruban d'étanchéité, échantillons et capuchons avec des septums préassemblés	1

Divers

Élément	Quantité dans l'expédition
 Kit d'évacuation, système Vanquish, avec : Pièce en L (coudée, paquet de 4, tailles différentes) Tube d'évacuation, longueur : 11,4 m Attaches (2 pièces) 	1
Kit de fusibles, système Vanquish Le kit comprend des fusibles 5AT pour la pompe, le passeur d'échantillon et le compartiment à colonnes (en option).	1
Schéma du flux	1
Échantillon de test cytochrome	1
Étiquette en papier pour le porte-étiquette magnétique	1
Tubes hélicoïdaux pour entourer les conduites ou les câbles	1
Porte-cartouche pour colonne de piégeage	1
Acclaim™ PepMap™ 100 C18 colonne HPLC (3 pièces)	1
Capuchon pour fermer les extrémités ouvertes des conduites de solvant (quand elles sont déconnectées de la pompe)	6
Bouchon pour fermer les ports de connexion ouverts dans la pompe (quand aucune conduite de solvant n'est connectée)	6

10.2.2 Compartiment à colonnes

Les pièces suivantes sont comprises dans le kit d'expédition du compartiment à colonnes.

Élément	Quantité dans l'expédition
Attaches pour colonne (paquet de 6)	1
Connecteurs d'ID de colonne (paquet de 2)	1
Câble System Interlink, longueur : 0,5 m Pour connecter le compartiment à colonnes au passeur d'échantillon.	1
Câble USB, de type A à type B, USB 2.0 à haut débit, longueur 1 m Pour connecter la pompe au compartiment à colonnes.	1

10.3 Accessoires en option

Boucles d'échantillonnage

Description	N° de référence
Boucle d'échantillonnage, biocompatible, MP35N, volume : 10 μL	6252.1960
Boucle d'échantillonnage, biocompatible, MP35N, volume : 25 μL (boucle d'échantillonnage standard)	6252.1940
Boucle d'échantillonnage, biocompatible, MP35N, volume : 100 μL	6252.1950

Portoirs d'échantillons et plaques à puits

Description	N° de référence
Portoir d'échantillons, accueillant 54 échantillons d'un D.E. de 12 mm	6850.1023
Avec code-barres 2D pour identification du type de portoir Vanquish	
Portoir d'échantillons, accueillant 96 échantillons d'un D.E. de 6 mm	6850.1026
Avec code-barres 2D pour identification du type de portoir Vanquish	
Portoir d'échantillons, accueillant 96 échantillons d'un D.E. de 7 mm	6850.1030
Avec code-barres 2D pour identification du type de portoir Vanquish	
Portoir d'échantillons, accueillant 96 échantillons d'un D.E. de 8 mm	6850.1034
Avec code-barres 2D pour identification du type de portoir Vanquish	
Portoir d'échantillons, accueillant 16 échantillons d'un D.E. de 15 mm	6851.1030
Avec code-barres 2D pour identification du type de portoir Vanquish	
Portoir d'échantillons, accueillant 9 échantillons d'un D.E. de 22,5 mm	6851.1020
Avec code-barres 2D pour identification du type de portoir Vanquish	
Pour obtenir des renseignements sur les commandes de consommables pour le passeur d'échantillon, tels que des plaques à puits, des tapis d'étanchéité, des échantillons et des bouchons, consultez thermofisher.com/ chromatographyconsumables.	

Vannes et accessoires du compartiment à colonnes

Élément	N° de référence
Connecteurs d'ID de colonne (paquet de 2)	6732.0610
Appareil de préchauffage passif, 0,1 mm x 680 mm, MP35N	6732.0184
Vanne 2 p-6 p, faible disp., 150 MPa, biocompatibilité	6250.1520
Vanne, 2 p-6 p, faible disp., VICI, C82	6041.0004B
Vanne, 2 p-6 p, faible disp., 1/16 po, 0,25 mm, PAEK	6821.5120
Vanne, 2 p-6 p, faible disp., 1/32 po, 0,1 mm, PAEK	6820.6232
Bouchon de raccordement pour vanne, 2 p-6 p, faible disp., 1/32 po, 0,1 mm, PAEK (paquet de 6)	6820.1320
Bague de verrouillage de vanne VICI	6253.1651

Isolateur USB

Élément	N° de référence
Isolateur USB L'isolateur USB améliore la fiabilité de la connexion USB entre le port USB d'un module du système HPLC et l'ordinateur du système de données.	6287.0540

Paillasses et stabilisateurs du système

Élément	N° de référence
Kit de serrage de paillasse Stabilise les piles du système avec une hauteur maximale de 139 cm	6036.1740
Kit de stabilisation du système empilé Stabilise les piles du système avec une hauteur maximale de 123 cm	6036.1710
lonBench avec kit de montage empilé Stabilise les piles du système avec une hauteur maximale de 123 cm	6036.1720

Flux de travail

Description	N° de référence
Kit de flux de travail Tandem pour injection directe, Vanquish Neo	6250.1030
fluidique du flux de travail d'injection directe Tandem Vanquish Neo. En plus du kit, le flux de travail requiert les modules du système suivants :	
 1 système Vanquish Neo (configuration standard) 	
 1 compartiment à colonnes Vanquish Neo avec 2 vannes à faible dispersion (2 à 6 p, 150 MPa) (n° de référence 6250.1520) 	
 1 pompe binaire Vanquish Neo 	
Pour obtenir des informations sur la configuration du flux de travail d'injection directe Tandem Vanquish Neo, consultez le <i>Guide de l'utilisateur du système Vanquish Neo</i> , révision 2.0 ou ultérieure.	

Divers

Élément	N° de référence
Kit d'application standard Vanquish Neo	6250.1010
Kit pour la réinstallation et la requalification d'un système Vanquish Neo utilisé. Il comprend :	
 Capillaire colonne – détecteur, nanoViper, gainage en silice fondue/PEEK, 1/32 po, D.I. x D.E. x longueur 20 μm x 280 μm, pour les applications nano LC 	
Capillaire colonne – détecteur, silice fondue/PEEK, D.I. x D.E. x longueur 50 μm x 280 μm, pour les applications capillaires et micro LC	
 Kit d'expédition de capillaire comprenant : 	
 Capillaire colonne – détecteur, silice fondue, D.I. x D.E. x longueur 20 μm x 280 μm x 1,5 m, pour les applications nano LC 	
 Capillaire colonne – détecteur, silice fondue, D.I. x D.E. x longueur 50 μm x 280 μm x 1,5 m, pour les applications capillaires et micro LC 	
 Capillaire, nanoViper, gainage en silice fondue/PEEK, 1/32 po, D.I. x D.E. x longueur 20 μm x 360 μm x 550 mm, pour les applications nano LC (2 pièces) 	
 Capillaire pompe – passeur d'échantillon, nanoViper, gainage en silice fondue/PEEK, 1/32 po, D.I. x D.E. x longueur 50 μm x 360 μm x 350 mm, pour les applications capillaires et micro LC 	
 Capillaire colonne – détecteur, nanoViper, gainage en silice fondue/PEEK, 1/32 po, D.I. x D.E. x longueur 50 μm x 360 μm x 550 mm, pour les applications capillaires et micro LC (4 pièces) 	
 Coupe-tube pour les tubes en silice fondue 	
• Acclaim™ PepMap™ 100 C18 colonne HPLC (3 pièces)	
Porte-cartouche pour colonne de piégeage	
 Échantillon de test cytochrome 	

10.4 Consommables et pièces de rechange

Pour obtenir des informations sur les numéros de référence des pièces de rechange spécifique à des flux de travail, consultez le *Guide de l'utilisateur du système Vanquish Neo,* révision 2.0 ou ultérieure.

10.4.1 Capillaires, accessoires des capillaires et évacuation

Capillaires

Description	N° de référence
Capillaire pompe – passeur d'échantillon, nanoViper, silice fondue, D.I. x longueur 20 μm x 350 mm	6250.5240
Capillaire passeur d'échantillon – compartiment à colonnes/ colonne, nanoViper, gainage en silice fondue/PEEK, 1/32 po, D.I. x longueur 20 μm x 550 mm	6250.5260
Capillaire pompe – passeur d'échantillon, nanoViper, gainage en silice fondue/PEEK, 1/32 po, D.I. x longueur 50 μm x 350 mm	6250.5540
Capillaire passeur d'échantillon – compartiment à colonnes/ colonne, nanoViper, gainage en silice fondue/PEEK, 1/32 po, D.I. x longueur 50 μm x 550 mm	6250.5560
Capillaire colonne – détecteur, nanoViper, gainage en silice fondue/PEEK, 1/32 po, D.I. x D.E. x longueur 20 μm x 280 μm x 1 m, pour les applications nano LC	6041.5292
Capillaire colonne – détecteur, nanoViper, gainage en silice fondue/PEEK, 1/32 po, D.I. x D.E. x longueur 20 μm x 360 μm x 1 m, pour les applications nano LC	6041.5293
Capillaire colonne – détecteur, nanoViper, gainage en silice fondue/PEEK, 1/32 po, D.I. x D.E. x longueur 50 μm x 280 μm x 1 m, pour les applications nano LC	6041.5294
Capillaire colonne – détecteur, nanoViper, gainage en silice fondue/PEEK, 1/32 po, D.I. x D.E. x longueur 50 μm x 360 μm x 1 m, pour les applications nano LC	6041.5295

Accessoires des capillaires

Description	N° de référence
Bouchon de raccordement, Viper, biocompatible	6040.2303
Raccord-union Viper	6040.2304

Outils capillaires

Description	N° de référence
Coupe-tube pour les tubes en silice fondue	6720.0016
Coupe-tube pour les tubes PEEK	6300.0401

Description	N° de référence
Tournevis dynamométrique, Viper, pièce de base avec adaptateur d'extension et embouts pour serrer les capillaires nanoViper et le siège de l'aiguille du passeur d'échantillon	6250.2110

Kit d'évacuation

Description	N° de référence
Kit d'évacuation, système Vanquish, avec :	6036.1120
 Pièce en L (coudée, paquet de 4, tailles différentes) 	
 Tube d'évacuation, longueur : 6,25 m 	
 Attaches de tubes d'évacuation (paquet de 2) 	

10.4.2 Pièces de rechange et outils du système

Pièces de rechange du système

Description	N° de référence
Bac à solvants (sans insertion de bouteille)	6036.1350
Insertion de bouteille pour bac à solvants	6036.1371
Dispositif d'affichage Vanquish	6036.1180
Vanquish System Controller	6036.1130-01

Pour les pièces de rechange du contrôleur du système Vanquish, reportez-vous à la section Pièce de rechange du contrôleur du système (> page 419).

Outils

Description	N° de référence
Outil de verrouillage du socle du système	6036.1160
Tournevis dynamométrique, Viper, pièce de base avec adaptateur d'extension et embouts pour serrer les capillaires nanoViper et le siège de l'aiguille du passeur d'échantillon	6250.2110
Kit d'outils, VN, incluant :	6250.2100
 Tournevis type Torx T10, en forme de L 	
 Tournevis type Torx T20, en forme de L 	
 Tournevis type Torx T30 	
 Clé hexagonale, taille 6 	
 Outil de manipulation des joints (pompe) 	
 Tournevis dynamométrique, Viper, pièce de base avec adaptateur d'extension et embouts pour serrer les connecteurs nanoViper et le siège de l'aiguille du passeur d'échantillon 	
Coupe-tube pour les capillaires PEEK	
 Coupe-tube pour les capillaires en verre de silice fondue 	

10.4.3 Pièce de rechange du contrôleur du système

Description	N° de référence
Carte SD	6439.0034
Câble System Interlink, longueur : 0,5 m Pour connecter la pompe au passeur d'échantillon.	6310.2012
Câble System Interlink, longueur : 0,25 m Pour connecter les éléments suivants : • Socle du système au contrôleur du système • Contrôleur du système à la pompe	6310.2013

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

10.4.4 Solvants et systèmes de rinçage

Réservoirs et éléments des réservoirs

Description	N° de référence
Réservoir, 1 l, avec bouchon	2270.0012
Réservoir, 0,25 l, avec bouchon	2270.0026
Bouchon pour réservoirs, bouchon à visser (paquet de 4)	6270.0013
Capuchon protecteur pour refermer les trous ouverts dans le bouchon de réservoir (paquet de 20)	6000.0047
 Guide de maintien (diamètre de 1,6 mm) pour garder les conduites suivantes dans le bouchon de réservoir (paquet de 5) : Conduites de solvant (entre le réservoir et la pompe) Conduites de rinçage (entre le réservoir et le dégazeur dans la pompe et entre le réservoir et la pompe de rinçage du passeur d'échantillon) 	6261.0216

Description	N° de référence
Guide de maintien (diamètre de 3 mm) pour garder la conduite de rinçage des joints dans le bouchon de réservoir (paquet de 5)	6000.0042
 Kit de capuchons et guides de maintien (diamètre de 1,6 mm) pour bouchons de réservoir de liquide de rinçage des joints, avec : Capuchon protecteur pour refermer les trous ouverts dans le bouchon de réservoir (paquet de 10) 	6030.9101
 Guide de maintien pour conduite de réservoir de liquide de rinçage des joints pour garder la conduite de rinçage des joints en place dans le bouchon du réservoir (paquet de 5) 	

Conduites de solvant, conduites de réservoirs de rinçage et filtres de conduites de solvant

Description	N° de référence
Kit de solvant et de conduites de rinçage pour le rinçage intérieur et extérieur de l'aiguille, y compris des conduites de réservoirs et étiquettes pour solvants	6250.2305
Conduites de solvant (A, B) entre les réservoirs de solvant et l'orifice d'admission de la pompe	6251.0096
Le kit de tubes à faible pression pour la pompe comprend aussi des conduites de solvant. Reportez-vous à la section Pièce de rechange de la pompe (> page 421).	
Filtre de conduite de solvant, porte-filtre (paquet de 6) (frittés de filtre ou adaptateur de conduite de solvant non compris)	6268.0115
Adaptateur de conduite de solvant (1/8 – 1/16) (paquet de 6)	6850.1056
Filtre de conduite de solvant, fritté de filtre, biocompatible, 10 μm (paquet de 10)	6268.0117

Tubes et éléments des dispositifs de rinçage

Description	N° de référence
Kit de tubes péristaltique et de rinçage	6044.1150
Ce kit inclut un tube péristaltique (PharMed), une tubulure en silicone et des raccords de tube à utiliser pour le	
 Système de rinçage des joints dans la pompe et dans le passeur d'échantillon 	
 Pompe d'évacuation dans le passeur d'échantillon 	
AVIS : Utilisez la tubulure en silicone épaisse (et raccords de tube correspondants) dans le passeur d'échantillon. Utilisez la tubulure en silicone mince (et raccords de tube correspondants) dans la pompe. Utilisez les raccords à visser pour la tête de pompe.	
Détecteur de rinçage des joints (détecteur de gouttes) avec pied	6044.1898A

Description	N° de référence
Kit de vis et d'étanchéité arrière VN (pompe VN), plaque de rinçage des joints/tête de pompe, incluant :	6251.2234
 Joint de la tête de pompe, pour la plaque de rinçage des joints (joints toriques 11x1,5) (2 joints) 	
 Joint de la tête de pompe, pour la vis de la plaque de rinçage des joints (joints toriques 6,5x1) (8 joints) 	
 Vis, plaque de rinçage des joints (8 vis) 	
Joint de piston (RP) et joint de rinçage des joints, tête de pompe (paquet de 2 pour chacun)	6251.1143
Kit PEEK de rinçage de l'aiguille, incluant	6252.1930
 Capillaire PEEK d'évacuation de l'orifice de rinçage (2 pièces) 	
 Connexion entre l'orifice de refoulement de la pompe de rinçage et l'orifice de rinçage 	

10.4.5 Pièce de rechange de la pompe

Tête de pompe et composants

Description	N° de référence
Tête de pompe (bloc complet)	6251.1102
Piston	6251.1155
Outil de manipulation des joints	6251.5122
Joint de piston (RP) et joint de rinçage des joints, tête de pompe (paquet de 2 pour chacun)	6251.1143
Bague de support (paquet de 2)	6040.0012
Clapet de non-retour, clapet de non-retour d'aspiration	6251.1153
Clapet de non-retour, clapet de non-retour de refoulement	6044.2310
Kit de capillaires, tête de pompe, incluant : • Tube en U • Capillaire entre la tête de pompe et le débitmètre	6251.0086
 Kit de vis et d'étanchéité arrière VN (pompe VN), plaque de rinçage des joints/tête de pompe, incluant : Joint de la tête de pompe, pour la plaque de rinçage des joints (joints toriques 11x1,5) (2 joints) 	6251.2234
 Joint de la tête de pompe, pour la vis de la plaque de rinçage des joints (joints toriques 6,5x1) (8 joints) 	
 Vis, plaque de rinçage des joints (8 vis) 	

Débitmètre

Description	N° de référence
Kit de filtre, débitmètre, incluant vis pour filtre, joint pour vis et fritté de filtre (2 μm)	6251.0111
Pour remplacer le filtre ou le fritté de filtre dans les deux entrées du débitmètre, deux kits sont nécessaires.	
Conduite d'évacuation du débitmètre, reportez-vous à <i>Kit de tubes faible pression</i>	

Kit de tubes faible pression

Description	N° de référence
Kit de tubes faible pression, pompe, incluant : • Conduite d'évacuation du débitmètre entre le débitmètre et l'orifice de refoulement de l'évacuation	6251.0150
 Conduites de solvant (A, B) entre les vannes d'arrêt et les têtes de pompe Conduites de solvant entre le dégazeur et la vanne d'arrêt 	
Reportez-vous aussi à la section Solvants et systèmes de rinçage (> page 419) et Capillaires, accessoires des capillaires et évacuation (> page 417).	

Kit de maintenance

Description	N° de référence
Kit de maintenance de la pompe, incluant :	6251.0130
 Vis, plaque de rinçage des joints (8 vis) 	
 Tubes en silicone (tubes transparents) 	
 Tube péristaltique (PharMed, tube blanc) 	
 Raccord de tube (droit) 	
 Raccord de tube (angle 90°) 	
 Raccord de tube (D.I. 1/16 po) (paquet de 4) 	
 Joint de piston (paquet de 4) 	
 Joint de la tête de pompe, pour la plaque de rinçage des joints (joints toriques 11 x 1,5) (paquet de 2) 	
 Joint de la tête de pompe, pour les vis de la plaque de rinçage des joints (joint torique 6,5 x 1) (paquet de 8) 	
 Joint de rinçage des joints (paquet de 4) 	
 Fritté de filtre de conduite de solvant (biocompatible, 10 μm) (paquet de 10) 	

Divers

Description	N° de référence
Capillaire nanoViper, silice fondue, D.I. x longueur 20 μm x 350 mm, à utiliser comme capillaire de contre-pression pendant la maintenance de la pompe	6250.5240
Bouchon de raccordement, Viper, biocompatible	6040.2303
Kit de tubes et seringues, incluant : • Seringue en plastique de 12 ml (paquet de 5) • Tubes en silicone (D.E. x D.I. 2,8 x 1,3), 3 m	6000.0010
Tubes de rinçage, silicone (D.E. x D.I. 2,8 x 1,3), 3 m	6007.9100
Écouvillons de nettoyage (paquet de 25)	6040.0007
Kit de fusibles, système Vanquish Pour la pompe, le passeur d'échantillon et le compartiment à colonnes, utilisez des fusibles à fusion lente 5 AT, 230 V c.a.	6036.0002

10.4.6 Pièces de rechange du passeur d'échantillon

Capillaires

Description	N° de référence	
Bouchon de raccordement, Viper, biocompatible	6040.2303	
Capillaire nanoViper, D.I. x L 50 μm x 120 mm	6252.5124	
Capillaires d'évacuation avec pièce en T	6252.1930	
Pour des capillaires supplémentaires, reportez-vous à la section Capillaires, accessoires des capillaires et évacuation () page 417).		

Composants fluidiques

Description	N° de référence
Kit de capillaires pour les composants fluidiques internes du passeur d'échantillon, incluant	6252.1920
 Capillaire Viper, entrée du capteur de pression 	
Capillaire Viper, sortie du capteur de pression	
Capillaire Viper, dispositif de mesure	
Capillaire Viper, clapet de non-retour	
 Capillaire PEEK d'évacuation de l'orifice de rinçage (2 pièces) 	
 Capillaire nanoViper, blindé FS/PEEK, 1/32 po 	
Tête de dispositif de mesure, volume 100 μL	6252.1753
Siège de l'aiguille	6252.2470
Unité d'aiguille	6252.1130
Clapet de non-retour, incluant le capillaire	6252.2310

Description

N° de référence

Pour les tubes et les dispositifs de rinçage, reportez-vous à la section Solvants et systèmes de rinçage (> page 419).

Divers

Description	N° de référence
Kit de portes avant, comprenant une porte droite et une porte gauche	6850.0100
Couvercle isolant	6252.1647
Verrou d'expédition pour carrousel	6850.5849
Tournevis dynamométrique pour serrer les raccords nanoViper et le siège de l'aiguille	6250.2110
Tubes de rinçage, silicone (D.E. x D.I. 2,8 x 1,3), 3 m	6007.9100
Kit de fusibles, système Vanquish	6036.0002
Pour la pompe, le passeur d'échantillon et le compartiment à colonnes, utilisez des fusibles à fusion lente 5 AT, 230 V c.a.	

10.4.7 Pièce de rechange du compartiment à colonnes

Description	N° de référence
Attaches pour colonne (paquet de 6)	6722.0290
Couvercles (paquet de 2)	6253.1647
Kit de portes avant, comprenant une porte droite et une porte gauche	6250.1620
Joint de rotor, 2 p-6 p, C72/C82, vanne E5 1/16 po, 0,15 mm	6041.0008B
Joint de rotor, 2 p-6 p, vanne C2/C3, 1/16 po, 0,25 mm	6820.0035
Stator, 2 p-6 p, vanne C72/C82 1/16 po 0,15 mm N60C	6041.0007
Stator, 2 p-6 p, vanne C1/C2 1/16 po 0,25 mm	6820.0034
Kit de fusibles, système Vanquish Pour la pompe, le passeur d'échantillon et le compartiment à colonnes, utilisez des fusibles à fusion lente 5 AT, 230 V c.a.	6036.0002

10.4.8 Câbles d'interface et cordons d'alimentation

Câbles d'interface

Description	N° de référence
Câble System Interlink, longueur : 0,5 m Pour connecter la pompe au passeur d'échantillon.	6310.2012
Câble System Interlink, longueur : 0,25 m Pour connecter les éléments suivants : • Socle du système au contrôleur du système • Contrôleur du système à la pompe	6310.2013
Câble USB, de type A à type B, USB 2.0 à haut débit, longueur : 0,5 m Pour connecter la pompe au passeur d'échantillon.	6720.8910A
Câble USB, de type A à type B, USB 2.0 à haut débit, longueur, 1 m Pour connecter la pompe au compartiment à colonnes.	6035.9035A
Câble USB, de type A à type B, USB 2.0 à haut débit, longueur, 5 m Pour connecter la pompe à l'ordinateur.	6911.0002A
Câble de signalement Digital I/O, 6 broches, longueur de câble : 5 m	6036.0006

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 et l'utilisation des ports Digital I/O (entrée/sortie numériques).

11.1 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.

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 :

E Su	e logo vert est apposé sur les dispositifs qui ne comportent pas les ubstances dangereuses répertoriées par la réglementation.
Le ap ré d' us er ris Pc w	e logo orange, qui encercle une valeur à un ou deux chiffres, est pposé sur les dispositifs qui comportent des substances dangereuses épertoriées par la réglementation. Le nombre indique la durée 'utilisation sans risques pour l'environnement (environment-friendly se period, EFUP) du dispositif. Pendant cette période, le dispositif (si mployé conformément à son utilisation prévue) ne présente pas de isques graves pour la santé humaine ou l'environnement. our 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 232: 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	
2.0	VN-A10-A-02, VN-C10-A-01, VN-P10-A-01, VN-S10-A-01	
1.0	VN-A10-A-02, VN-C10-A-01, VN-P10-A-01, VN-S10-A-01	

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

11.2 E/S numériques (affectation des broches)

Les ports E/S numériques (Dig I/O) peuvent être utilisés pour échanger des signaux numériques avec des appareils externes, par exemple, un spectromètre de masse. Chaque port comprend :

- une entrée numérique
- une sortie relais



Illustration 233: Port d'entrée et sortie numériques

Broche	Description – Nom du signal		
1	Non utilisé		
2	Sortie relais — Relay_NC (contact normalement fermé)		
3	Masse — GND		
4	Entrée numérique — Input		
5	Sortie relais — Relay_COM COM est le contact habituel pour NO et NC. Si le relais n'est pas activé ou bien si l'appareil est hors tension, la connexion s'effectue entre COM et NC. Si le relais est activé, la connexion s'effectue entre COM et NO.		
6	Sortie relais — Relay_NO (contact normalement ouvert)		

Broche	Couleur du fil	Désignation du signal	Niveau du signal	Remarques
1	Rose			Non utilisé
2	Gris	Sortie relais — Relay_NC	Contact sec 0-24 V, 0-100 mA	Contact d'ouverture
3	Verte	Masse — GND	Masse	Potentiel de référence
4	Jaune	Entrée numérique — I nput	Entrée (faible activité) : Activée : 0-0,4 V Désactivée : 2,2-5 V Résistance de rappel vers le haut : 47 kΩ à 5 V	 Entrée numérique ; le potentiel de référence est la masse. Veuillez noter les points suivants : La tension en entrée maximale ne doit pas dépasser +5 V par rapport à la masse. La tension en entrée minimale ne doit pas être inférieure au potentiel de la masse.
5	Blanc	Sortie relais — Relay_COM	Contact sec	Contact habituel de NO et NC
6	Marron	Sortie relais — Relay_NO	Contact sec 0-24 V, 0-100 mA	Contact de fermeture

Le tableau suivant répertorie les fonctions affectées aux broches du connecteur ainsi que la couleur du fil branché sur chaque broche.

Consultez également

Connexion d'un spectromètre de masse () page 118)
Index

Α

additifs	
compatibilité générale	144
utilisation	145
algues	145
alimentation	
boutons	62, 136
commutateurs	62, 136
exigences	398
mettre en marche	136, 148
alimentation électrique	79 <i>,</i> 80
appareil de préchauffage passif	
installation	92
remplacement	313
appareil LAN	
intégration du système	138
arrêt	
courte durée	177
longue durée	177
arrêt de courte durée	177
arrêt de longue durée	177
-	

В

bac à solvants	56
insertion de l'insertion de bouteille	330
remplacement	326
retrait de l'insertion de bouteille	329
bac d'évacuation	126
bague de support	
installation	206
remplacement	204
retrait	204
balayage d'inventaire	147
Barre à LED 149, 153,	354
bouchons de flacons	167
bouton ROTATE	165
boutons du panneau de commande	150

С

câble LAN	
raccordement	115
câbles System Interlink	
raccordement	113
câbles USB	

raccordement	110
capillaires	
guide	120
installation	121
capillaires nanoViper	
connecter	122
caractéristiques	
dimensions	399
performance de la pompe	402
performance du compartiment à colonne	<u></u>
performance du compartiment à colonne.	, 405
caractéristiques techniques	207
avigances relatives à l'alimentation électri	iquo
exigences relatives a railmentation electric	ique
	398
niveau de pression acoustique	399
performance du contrôleur du système	404
performance du passeur d'échantillon	403
performance du système	401
poids	400
carrousel 49) , 50
chargement	165
code couleurs	. 50
faire pivoter	165
porte-flacon supplémentaire	. 50
carte SD	
remplacement	321
changement	
liquide de rincage de l'aiguille	172
liquide de rincage des joints	172
changement de type de solvant 163	172
clanet de non-retour	
rincage de l'aiguille	281
tâte de nomne	201
clanat de pompe	220
nottovago	220
nettoyage	229
remplacement du clapet d'aspiration	220
remplacement du clapet de refoulement	228
clapet de non-retour d'aspiration	
nettoyage	229
remplacement	226
clapet de non-retour de refoulement	
nettoyage	229
remplacement	228
clapet de non-retour de rinçage de l'aiguille	
remplacement	281

classe de sécurité	26
cloison	167
code-barres (Vanquish)	51, 167
colonnes	
installation	94
remplacement	303
compartiment à colonnes	
caractéristiques	35
composants internes	53
installation	86
préparation du retrait	341
principe de fonctionnement	54
retrait du module à glissières	342
thermostat	175
compartiment à échantillons	49
carrousel	50
thermostat	50
composants internes	175
compartiment à colonnes	E 2
passour d'échaptillen	
	44 11
pompe	41
concentration en chiorure	145
condensation	16, 320
conditions d'entreposage	
humidite	398
température	398
conditions d'utilisation	81
altitude	398
chocs mécaniques	82
humidité	81, 398
interférence électromagnétique	82
température	81, 398
ventilation	82
vibrations	82
conduites d'évacuation	124
conduites de solvant	
entre la vanne d'arrêt et la tête de pon	npe
	242
entre le réservoir et la pompe	237
maintenance	237
vide	237
Conformité aux normes UL/CSA	429
conformité avec la réglementation	
connecteurs	
contrôleur du système 1	07 318
modules du système	106
connecteurs des modules du système	106 106
connecteurs du contrôleur du système 1	100 07 210
connecteurs au controleur au systeme. 1	57, 510

connexions à l'alimentation électrique consignes	115
installation	76
maintenance	181
utilisation	142
consignes de sécurité	172
cas d'urgence	31
cordon d'alimentation	91 80
entretien	181
équipement de protection	101
généralités	27 26
installation	20
maintenance	121
qualification du personnel	101
	····· 27 20
sécurité électrique	29 20
utilisation	1/2
contrôleur du système	142 //
connector 10	40
data	155
houro	155
installation	133
maintananco	515
	314 155
romplacement de la carte SD	100
retrait	521 214
retrait	314
version du microprogramme	155
cordon d'alimentation	0, 320
	97,98
installation	00
IIIstallation	98
retrait	97

D

déballage du compartiment à colonnes	. 72
déballage du système	. 70
débit (interdépendance avec la pression de la	1
pompe)	169
débitmètre	
maintenance	232
remplacement de la conduite d'évacuation	n
	234
remplacement des filtres	232
décontamination	191
DEEE	429
dépannage	353
alarme de fuite	395
analyse instable	385

faible reproductibilité3	85
généralités 3	54
mauvaise précision de l'aire des pics 3	84
messages	56
pics fantômes 3	83
pics négatifs 3	83
pics parasites	84
problèmes d'alimentation 3	92
problèmes de communication	93
problèmes de forme de pics	82
problèmes de la ligne de base 3	89
problèmes de la régulation de la températu	re
du compartiment à colonnes 3	91
problèmes de la régulation de la températu	re
du passeur d'échantillon 3	91
problèmes de pics 381, 3	83
problèmes de pression	86
scripts	79
temps de chargement prolongé	90
variation des temps de rétention	88
dépannage du système	79
détecteur de gouttes (remplacement)	98
détecteur de rinçage des joints (remplacement)
	, 98
détection de fuites	57
déverrouillage du socle du système	73
Dig I/O 4	31
dimensions	99
dispositif d'affichage	
déplacement de la position frontale à la	
position latérale 1	60
désinstallation	23
installation	99
maintenance 3	14
positions	59
raccordement1	11
réglage1	59
remplacement	23
dispositif de rincage de l'aiguille	
conduite de sortie de la pompe de rincage	
	75
conduites de réservoir 273. 2	77
purge	33
remplacement des conduites 273-275-2	77
remplissage des réservoirs	31
rincage de l'aiguille	74
donnees techniques	
altitude de fonctionnement	98

dimensions	399
exigences relatives à l'alimentation électri	que
	398
humidité d'entreposage	398
humidité de fonctionnement	398
niveau de pression acoustique	399
poids	400
température d'entreposage	398
température de fonctionnement	398

Ε

F

FCC 4	29
filtre de conduite de solvant	
assemblage 129, 2	44
maintenance 2	37
remplacement des filtres 2	43
remplacement du fritté 2	43
flacon	
rinçage de l'aiguille 1	31
rinçage des joints 1	34
solvant 1	29
fonction de vanne de captage 91, 3	10
fuite 3	95
fuites de liquide 3	95

fuseau horaire	157
fusibles	332
fusibles de l'alimentation principale	332

G

gaine de tubulure	120
gants	28
guides de tubulure	120

I

icônes dans l'interface utilisateur Vanquish
identification du type de portoir 51, 167
insertion de bouteille
insertion 330
retrait 329
installation
capillaires et tubes 120
conduites d'évacuation 124
connexions à l'alimentation électrique 115
consignes de sécurité
exigences relatives au lieu d'installation 78
système 83
Interface utilisateur Vanquish
fuseau horaire 157
licences d'utilisation 157
Option personnalisée 138
possibilité de connectivité 138
possibilité DHCP 138
réglages de connectivité 157
réglages de langue 157
Interlink
interrupteurs d'alimentation principaux 62, 136

J

joint de piston	
installation	212
remplacement	209
retrait	209
joint de rinçage des joints	
installation	218
remplacement	216
retrait	216
joint de tête de pompe	
installation	222
remplacement	220
retrait	220

joint du rotor de la vanne de commutation V	ICI
remplacement	311

Κ

kit d	'expédition	409
-------	-------------	-----

L

LED D'ÉTAT 14	19, 354
licences d'utilisation	157
liquide de fuite	124
liquide de rinçage de l'aiguille	131
changement	172
liquide de rinçage des joints	
changement	172
livraison	69
lunettes de sécurité	28

Μ

maintenance	179,	186
consignes de sécurité		181
contrôleur du système		314
décontamination		191
dispositif d'affichage		314
fusibles		332
introduction		180
nettoyage		191
portes		333
règles générales		184
tester l'étanchéité		395
maintenance annuelle		189
maintenance hebdomadaire		187
maintenance régulière		188
Marquage CE		428
Marquage cTUVus		429
Marquage EAC		428
Marquage RoHS		428
marquage UKCA		429
matériel compris		. 69
messages		
compartiment à colonnes		378
généralités		356
messages communs		356
passeur d'échantillon		373
pompe		362
volet notification		354
messages concernant la pompe		362
messages d'avertissement	•••••	. 22

mettre le système en marche 14	8
mettre le système sous tension	
mise en marche initiale 13	6
régulière 14	8
mise en marche initiale du système 13	6
mise sous et hors tension 6	52
mode d'injection avec division de flux 4	-5
module à glissières	
installation 34	5
retour 34	4
module à glissières (pompes)	
retrait	57

Ν

nettoyage	191
niveau de pression acoustique	399

0

0	otimiser	les	performances		169
	Summer	105	periormanees	••••••	105

Ρ

couvercles 1	67
orientation 1	66
plaque signalétique	24
plaques signalétiques	23
poids 4	00
poignées	
fixation au système 3	50
poignées de manutention	
fixation au système 3	50
pompe	
caractéristiques	34
composants internes	41
préparation du retrait 3	36
principe de fonctionnement	42
retrait du module à glissières	37
pompe (rinçage) 189, 1	90
porte	
décrochage 3	33
fermeture	85
mécanisme d'ouverture du passeur	
d'échantillon	85
mécanisme de fermeture du passeur	
d'échantillon	85
ouverture	85
remplacement 3	33
retirer 1	80
retrait1	80
portoir d'échantillons	50
code-barres	51
orientation 1	66
présentation (fonctionnelle)	33
pression de la pompe (interdépendance avec le	e
débit) 1	69
principe de fonctionnement	
compartiment à colonnes	54
passeur d'échantillon	45
pompe	42
purge	65

R

raccordement	
cordon d'alimentation	320
raccordements fluidiques	
conduites d'évacuation	124
dispositif d'évacuation	124
tubes de la pompe d'évacuation	268
Ready state	176
Ready Temp Delta	176

réglage de la position du dispositif d'affichage

réglage du décalage du capteur de débit 173
réglages de connectivité 157
réglages de langue 157
réinitialiser aux valeurs par défaut d'usine 380
réparation en atelier 185
réponse d'injection 118
réservoir
rinçage de l'aiguille 131
rinçage des joints 134
solvant 129
réservoirs de solvant 129, 146
retour
module à glissières 344
système 185
rinçage
après utilisation 146
pompe 189, 190
rinçage faible 131
rinçage fort 131
ruban d'étanchéité 167

S

script de réinitialisation des réglages d'usine pa	ar
défaut 3	380
Script Démarrage automatique 139, 1	61
Script Nettoyage du système 3	379
Script Tester la contre-pression du système 3	379
Script Tester le serrage du système 3	379
Script Initialiser réglage du système 3	380
serre-câbles 1	17
socle du système	38
déverrouillage	73
verrouillage 1	19
solvant	
changement 1	46
compatibilité générale 1	44
concentration en chlorure 1	L45
concentration tampon 1	L45
étalonnage 1	170
plage pH 1	44
qualité et filtration 1	L45
utilisation 1	L45
spectromètre de masse (connexion) 1	18
stator de la vanne de commutation VICI	
remplacement3	311
support de tubulure 1	120

symboles de sécurité 2	2, 23
System Interlink 39	, 106
système	
arrêt	. 177
caractéristiques	34
dépannage	. 379
description	37
équilibrage	. 163
état	. 155
optimiser les performances	. 169
préparation du transport	. 347
préparation pour l'utilisation	. 161
préparation pour l'utilisation initiale 139	, 161
vue d'ensemble	37
système de rinçage des joints 60), 282
conduite d'admission de la pompe de rin	çage
des joints	. 291
conduite d'admission de rinçage des join	ts
	. 288
conduite du réservoir de liquide de rinça	ge
des joints	. 284
conduites de rinçage des joints dans la	
pompe	. 296
détecteur de rinçage des joints	. 298
liquide de rinçage des joints	. 134
maintenance	. 282
pompe de rinçage des joints	. 293
remplissage de réservoir	. 134
rinçage	. 137
tester l'étanchéité	. 282
tube péristaltique	. 293
système de rinçage des joints arrière	60
système de rinçage des joints de piston	60

Т

tampons	
concentration	145
utilisation	145
tapis de couverture	167
Temps d'équilibrage	176
test d'étanchéité	
étanchéité du système	395
système de rinçage des joints	282
tubes de la pompe d'évacuation	268
test de fuite détaillé (pompe)	396
tête de pompe	
clapet de non-retour 226,	229
clapet de non-retour d'aspiration	226

clapet de non-retour de refoulement 228
installation 198
maintenance 195
remplacement 197
retrait 197
vue d'ensemble des éléments 196
tête du dispositif de mesure
connexion des capillaires
déconnexion des capillaires
remplacement
thermostatisation du compartiment à
échantillons 50
tournevis dynamométrique 122
transport
préparation du système 347
trou de guidage 120
tube péristaltique
remplacement 294
tubes 121
tubes de la pompe d'évacuation 268
recherche d'obstruction
remplacement 270
test
tester l'étanchéité 269
type de solvant
changement 163, 172

U

unité d'aiguille	
déverrouillage 25	52
position d'entretien 25	51
verrouillage 25	51
Universal Serial Bus 10)6
USB 10)6
utilisation 14	1
consignes de sécurité 14	2
consignes générales 14	7
éléments de commande 14	9
Interface utilisateur Vanquish	54
interruption pour une courte durée 17	7
interruption pour une longue durée 17	7
système de gestion de données	
chromatographiques6	54
utilisation prévue 2	25

vanne d'injection attribution des orifices 48 vanne de commutation pour colonne 54 fuite 304 vanne de commutation Vanquish..... 55 installation 88 remplacement 305 Vanne de commutation VICI...... 55 installation 90 remplacement 308 vanne de solvant attribution des orifices 49 vannes du passeur d'échantillon attribution des orifices 48 Vanquish User Interface icônes 155, 156 scripts 379 verrouillage du socle du système 39 vêtements de protection 28 vis de verrouillage d'expédition serrage...... 251 volet notification 354 voyant d'état Barre à LED 149, 354

V

valeurs par défaut d'usine

www.thermofisher.com

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

