Nicolet Summit LITE

Spectromètre FTIR



MANUEL DE L'UTILISATEUR

269-345500 Révision C Avril 2021

© 2020 Thermo Fisher Scientific Inc. Tous droits réservés.

Pour obtenir une assistance technique, contactez : www.thermofisher.com.

Thermo Fisher Scientific Inc. fournit cette documentation à ses clients à l'achat d'un produit. Cette documentation est protégée par copyright et toute reproduction intégrale ou partielle de celle-ci est formellement interdite, sauf autorisation écrite de Thermo Fisher Scientific Inc.

Le contenu de cette documentation peut faire l'objet de modifications sans préavis. Toutes les informations techniques sont fournies à titre de référence uniquement. Les configurations et spécifications du système du présent document remplacent toutes les informations précédentes.

Thermo Fisher Scientific, Inc. n'affirme en aucun cas que cette documentation est complète, précise ou exempte d'erreurs, décline toute responsabilité et ne peut être tenue pour responsable de toute erreur, omission, perte ou dommage causés par l'utilisation de la présente documentation, même si les consignes qu'elle contient sont suivies scrupuleusement.

Ce document ne fait pas partie du contrat de vente entre Thermo Fisher Scientific Inc. et l'acheteur. Ce document ne régit pas ni ne modifie de quelque manière les Conditions de vente, lesquelles régissent la résolution de tous les conflits pouvant survenir entre ces deux documents.

Usage exclusivement réservé à la recherche. Cet instrument ou accessoire n'est pas un dispositif médical et n'est pas conçu pour être utilisé pour la prévention, le diagnostic, le traitement ou la guérison de maladies.

AVERTISSEMENT



Évitez tout risque d'explosion ou d'incendie.

Cet instrument ou accessoire n'est pas conçu pour être utilisé dans une atmosphère explosive.

Table des matières

Bienvenue	1
Conventions employées	2
Commande de pièces	3
Nous contacter	3

Première utilisation du spectromètre Nicolet Summit LITE	5
Avant la livraison du spectromètre Summit	6
Sécurité personnelle dans le cadre de l'utilisation du spectromètre Summit	9
Déballage et première utilisation du spectromètre	16
Installation d'un accessoire d'échantillonnage	21
Comprendre la barre lumineuse du spectromètre	27
Consultation de vos données via l'application OMNIC Anywhere	30

Options et accessoires	31
Accessoires d'échantillonnage pour le spectromètre Summit	
Installation et entretien d'un kit de purge	

Maintenance	45
Calendrier de maintenance du spectromètre Summit	
Alignement des composants optiques du spectromètre	50
Calibration de la fréquence du laser	51
Nettoyage du spectromètre	52
Remplacement du desséchant	53
Remplacement de la source	59
Remplacement des fenêtres du compartiment à échantillons	63

Chapitre 1

Bienvenue

Le spectromètre infrarouge à transformée de Fourier (FTIR) Thermo Scientific Nicolet Summit LITE vous permet d'effectuer des analyses chimiques d'échantillons de substances en recueillant des données dans le domaine spectral de l'infrarouge moyen à l'aide de divers accessoires. Le système possède des fonctions de vérification intégrées, une puissante suite logicielle et de nombreuses autres caractéristiques qui facilitent l'acquisition des données. Vous pouvez installer du matériel en option et effectuer vous-même plusieurs procédures d'entretien et de maintenance. Le présent document, ou toute autre documentation fournie, contient les informations dont vous avez besoin.

Assurez-vous de lire le guide de sécurité fourni avec le système avant d'utiliser l'instrument.

Note Nous recommandons de constamment garder le spectromètre scellé, dessiqué et/ou purgé. Les dommages causés à l'équipement par un défaut d'étanchéité et de desséchant et/ou de purge ne sont pas couverts par la garantie. Si vous avez des questions concernant cette exigence, contactez-nous.

Conventions employées

Les mesures de sécurité et les autres informations importantes utilisent le format suivant :



Évitez tout danger. Indique une situation dangereuse qui, si elle n'est pas évitée, entraînera la mort ou des blessures graves.

AVERTISSEMENT



Évitez tout danger. Indique une situation dangereuse qui, si elle n'est pas évitée, peut entraîner la mort ou des blessures graves.

ATTENTION



Évitez tout danger. Indique une situation dangereuse qui, si elle n'est pas évitée, peut entraîner des blessures légères ou modérées.

À NOTER

Suivez les instructions précédées de cette mention afin d'éviter d'endommager le matériel système ou de perdre des données.

Note Contient des informations supplémentaires utiles.

Bienvenue

Commande de pièces

Pour commander des pièces, contactez-nous.

Si vous devez nous expédier l'instrument ou un accessoire pour réparation, appelez-nous ou envoyeznous un courriel au préalable pour connaître toute exigence relative à l'expédition, ou obtenir d'autres instructions.

Nous contacter

Pour obtenir une assistance technique, contactez : www.thermofisher.com.

[Cette page est intentionnellement laissée vide]

Première utilisation du spectromètre Nicolet Summit LITE

Pour vous familiariser avec votre nouveau spectromètre, le Nicolet Summit LITE, passez en revue les exigences relatives à la sécurité et au site d'installation de l'instrument avant l'arrivée de ce dernier. Le jour de la livraison, vous pourrez déballer et installer l'instrument ainsi que ses accessoires.

Avant la livraison du spectromètre Summit	6
Sécurité personnelle dans le cadre de l'utilisation du spectromètre Summit	9
Déballage et première utilisation du spectromètre	16
Installation d'un accessoire d'échantillonnage	21
Comprendre la barre lumineuse du spectromètre	27
Consultation de vos données via l'application OMNIC Anywhere	

Avant la livraison du spectromètre Summit

Préparez un espace de travail approprié pour le spectromètre avant livraison afin de garantir la précision des données et les performances à long terme de l'instrument.

Cette rubrique présente les facteurs environnementaux et électriques qui peuvent avoir une incidence sur les performances de l'instrument, et explique les actions requises à la livraison du spectromètre. Des explications plus complètes sur ces exigences sont disponibles à la rubrique « Informations relatives à la sécurité et au site d'installation » relative au spectromètre.

Déballage du spectromètre

Avant d'ouvrir la boîte, il convient de respecter deux étapes importantes à l'arrivée de l'instrument :

• Vérifiez si l'extérieur de la boîte d'expédition n'a subi aucun dommage

Si vous remarquez des signes de dommages, contactez notre service d'assistance technique ou votre distributeur local pour obtenir des instructions.

· Laissez le spectromètre atteindre la température ambiante

Dans la boîte d'expédition, le spectromètre est scellé dans un sac en plastique destiné à le garder au sec. Attendez 24 heures pour ouvrir le sac - le temps que le spectromètre atteigne la température ambiante. Si vous ouvrez le sac avant que le spectromètre ne se soit réchauffé, de la condensation pourrait se former dans l'instrument, au risque d'endommager ses composants optiques internes et de l'endommager de manière permanente.

La garantie ne couvre pas les dommages résultant de techniques de déplacement de l'instrument inappropriées ou du retrait de l'instrument hors du sac en plastique scellé avant que l'instrument n'ait atteint la température ambiante.

Préparation de l'espace de travail

Avant la livraison de l'instrument, veillez à ce que votre espace de travail réponde pleinement aux exigences relatives au site d'installation du spectromètre. En plus de prévoir un espace suffisant pour le spectromètre, prenez en compte plusieurs exigences environnementales et électriques.

Dimensions du spectromètre

Le spectromètre Summit ne requiert qu'un espace assez restreint, mais assurez-vous de laisser de la place tout autour de l'instrument pour que la chaleur qu'il génère puisse se dissiper par les aérations, et

pour vous permettre d'accéder facilement aux différents ports, au bouton d'alimentation et aux câbles de l'instrument.

- Poids de l'instrument Summit : 9,6 kg (21 lb)
- Dimensions (L x H x D) : 34 x 24 x 32 cm ; (13,3 x 9,6 x 12,7 po)

Facteurs environnementaux

Le spectromètre Summit est un instrument robuste, conçu pour être utilisé dans de nombreux environnements. Cependant, pour obtenir de meilleures performances, il convient de le conserver dans un environnement relativement exempt de poussière et d'humidité. Le spectromètre peut fonctionner de manière fiable entre 15 et 35 °C, mais ses performances sont optimales dans un environnement entre 20 et 22 °C.

L'humidité peut former une condensation à l'intérieur de l'instrument, susceptible d'endommager ses composants internes. Les quelques précautions suivantes sont généralement suffisantes pour protéger votre instrument de l'humidité :

- Veillez à ce que du desséchant soit toujours présent dans l'instrument, y compris pendant son stockage.
- Évitez les brusques changements de température.
 - Tenez l'instrument à l'écart des sources d'air froid ou chaud par exemple, des bouches de chauffage, des sorties de climatisation, ou des grandes fenêtres.

Si le spectromètre doit être conservé dans un environnement particulièrement humide, envisagez d'installer un kit de purge.

Conditions requises en matière d'alimentation électrique

L'alimentation électrique du spectromètre doit provenir de sources dédiées, ininterrompues, et exemptes des problèmes suivants :

- Coupures de tension
- Surtensions transitoires
- Changements de fréquence
- Autres perturbations de la ligne électrique.

Si vous suspectez des problèmes liés à l'alimentation électrique, nous vous recommandons d'effectuer un audit de la qualité de l'alimentation électrique du bâtiment. Pour plus d'informations, contactez notre service d'assistance technique ou votre fournisseur d'électricité.

Spécifications de l'alimentation électrique

Le tableau suivant énumère les spécifications requises en matière d'alimentation électrique. Pour toute question relative à ces spécifications, contactez le représentant du service après-vente de votre région.

Conditions requises	Spécifications
Courant d'entrée	1,5 A max.
Tension d'entrée	100 à 240 V CA
Fréquence du secteur	50 à 60
Perturbations de ligne	Les chutes de tension, les surtensions et les autres perturbations de ligne ne doivent pas dépasser 10 % de la tension d'entrée
Bruit	Moins de 2 V (mode commun) moins de 20 V (mode normal)

Sécurité personnelle dans le cadre de l'utilisation du spectromètre Summit

Bien que la conception du spectromètre Nicolet Summit LITE garantisse la sécurité d'exploitation de l'instrument, vous devez prendre certaines précautions pour vous protéger de dangers qui peuvent survenir lors de l'utilisation normale et de la maintenance de l'instrument.

ATTENTION



Sans être exhaustif, ce guide est une introduction aux dangers potentiels dont vous devez être conscient(e). Avant d'utiliser l'instrument, consultez la rubrique « Informations relatives à la sécurité et au site d'installation » pour obtenir une description complète des dangers potentiels.

Dangers potentiels associés à l'utilisation normale de l'instrument

Dans le cadre d'une utilisation normale de l'instrument, la plupart des dangers proviennent des sources suivantes :

- Échantillons et solvants potentiellement dangereux
- Exposition au séléniure de zinc (ZnSe)
- Parties du spectromètre chaudes ou qui génèrent de la chaleur
- Laser de l'instrument

Vous pouvez éviter de vous blesser ou d'endommager l'instrument en comprenant les dangers potentiels et en prenant certaines précautions.

Échantillons et solvants dangereux

Prenez des précautions particulières si vous prévoyez de mesurer des échantillons ou d'utiliser des solvants potentiellement dangereux, tels que des gaz sous pression ou des solvants corrosifs ou inflammables.

UTILISER UNE VENTILATION ADÉQUATE

Votre spectromètre ne nécessite pas de ventilation particulière, mais une ventilation supplémentaire peut être nécessaire avec certains types d'analyse. Assurez la ventilation adéquate de la salle de travail

si vous devez analyser des échantillons hautement toxiques, dissoudre des échantillons dans des solvants susceptibles d'interagir avec la source infrarouge, ou échantillonner des gaz inflammables.

La pyrolyse des solvants contenant des hydrocarbures halogénés peut produire de l'acide chlorhydrique (HCI), de l'acide fluorhydrique (HF) ou du phosgène (COCI₂).

AVERTISSEMENT



Évitez toute inhalation toxique. L'acide chlorhydrique, l'acide fluorhydrique et le phosgène sont des substances hautement toxiques. Si vous utilisez des solvants contenant des hydrocarbures halogénés, assurez-vous que votre zone de travail est correctement ventilée.

SOLVANTS VOLATILS ET INFLAMMABLES

La source infrarouge à l'intérieur du spectromètre peut enflammer des échantillons et des solvants inflammables et volatils. Prenez les mesures suivantes lorsque vous travaillez avec des échantillons et des solvants inflammables :

- Ne travaillez qu'avec les fenêtres d'isolation du compartiment à échantillons installées.
- Assurez-vous que l'espace de travail est correctement ventilé à l'aide d'un système de ventilation actif qui ne génère pas d'étincelles ou d'autres sources d'inflammation, et qui prévient l'accumulation de vapeurs inflammables dans l'atmosphère entourant l'instrument.
- Ne laissez pas de solvants ni d'échantillons inflammables à proximité de l'instrument.
- Ne laissez pas de solvants ni d'échantillons inflammables dans le compartiment à échantillons plus longtemps que nécessaire.
- Purgez le spectromètre avec de l'air propre et sec ou de l'azote.

SOLVANTS CORROSIFS

L'utilisation de solvants pouvant produire des vapeurs de HCI ou HF dans le compartiment à échantillons peut gravement endommager le système. Si vous utilisez des solvants halogénés, purgez l'instrument avec de l'air ou de l'azote propre et sec.

À NOTER

Votre garantie ne couvre pas les dommages causés à l'équipement en raison d'un défaut de purge de l'instrument.

Les vapeurs de HCI et de HF peuvent également compromettre le revêtement des fenêtres d'isolation du compartiment à échantillons en KBr. Si vous prévoyez de travailler régulièrement avec des solvants corrosifs, envisagez plutôt d'installer des fenêtres de compartiment à échantillons en ZnSe.

MATIÈRES RADIOACTIVES OU PRÉSENTANT UN DANGER BIOLOGIQUE ET AGENTS INFECTIEUX

Les échantillons biologiques tels que les tissus, les liquides corporels, les agents infectieux et le sang sont susceptibles de transmettre des maladies infectieuses. Suivez les protocoles du programme de biosécurité de votre organisation lorsque vous travaillez avec des matériaux potentiellement infectieux.

Matériaux toxiques

Le spectromètre Nicolet Summit OA comprend un cristal ATR en séléniure de zinc (ZnSe), et d'autres instruments Summit peuvent utiliser les fenêtres de compartiment à échantillons en ZnSe en option.

AVERTISSEMENT



Évitez toute inhalation et ingestion toxique.

Le séléniure de zinc (ZnSe) est toxique. Reportez-vous à la Fiche de données de sécurité du ZnSe du fabricant sur www.specac.com pour la manipulation et les contrôles d'exposition.

Sources de chaleur

Certaines parties du spectromètre peuvent devenir très chaudes dans le cadre d'une utilisation normale. Faites attention à la source infrarouge du spectromètre et aux aérations de l'instrument.



Figure 2-1 : Prudence à proximité de la source infrarouge et des aérations de l'instrument

La surface de la source infrarouge orientée vers l'extérieur est située sous l'instrument et peut devenir très chaude. Ne touchez pas la base de l'instrument lorsque ce dernier est en marche, ou à l'issue de son utilisation.

L'aération du spectromètre est effectué sur le côté gauche de l'instrument. Dans le cadre d'une utilisation normale, le spectromètre dissipe l'air chaud des composants internes vers la zone située à l'extérieur du spectromètre par ces trous d'aération. Laissez suffisamment d'espace libre autour des trous d'aération de l'instrument pour que l'air chaud puisse se dissiper.

Sécurité des lasers et des composants optiques

Vous ne serez jamais exposé(e) à des niveaux de rayonnement laser dangereux dans le cadre de l'utilisation normale du spectromètre. Cependant, si vous retirez le couvercle au cours d'une tâche d'entretien, vous devrez peut-être prendre des précautions particulières - par exemple, vous doter de lunettes de protection. Votre technicien d'entretien vous informera si cela est nécessaire.

AVERTISSEMENT



Évitez toute lésion corporelle.

Ne regardez jamais le faisceau laser ni sa réflexion. Ne manipulez ou ne modifier jamais le laser, même lors du remplacement d'un laser défectueux, sous peine de vous exposer à la lumière du laser ou à une tension élevée.

Dangers potentiels pendant la maintenance

La maintenance de l'instrument peut vous exposer à des risques autres que ceux rencontrés lors d'une utilisation normale. Au cours de la maintenance, la purge de l'instrument et le maniement de ses composants internes sont les principales sources de dangers potentiels.

Purge de l'instrument

Dans les environnements particulièrement humides, il est recommandé d'installer une source d'air propre et sec ou d'azote pour purger le spectromètre. La purge de l'instrument peut aider à protéger ses composants optiques internes des dommages provoqués par un environnement humide ou des solvants corrosifs, et garantir des résultats plus précis.

Des informations sur l'achat et l'installation d'un kit de purge pour le spectromètre sont disponibles à la rubrique <u>"Installation et entretien d'un kit de purge</u>".

DANGER Évitez tout risque d'incendie et d'explosion. Utilisez uniquement de l'azote ou de l'air sec pour purger le spectromètre. N'utilisez jamais de gaz inflammable, combustible ou toxique pour purger cet instrument. Le gaz de purge ne doit pas contenir d'huile ni d'autres matières réactives. La chaleur provenant de la source ou de l'absorption du laser peut enflammer des gaz inflammables ou des matières réactives dans le gaz de purge.

Manipulation de composants internes

En règle générale, il n'existe aucune raison justifiant le retrait du couvercle de l'instrument ou le remplacement des composants internes. Cependant, si vous devez retirer le couvercle dans le cadre de la maintenance, sachez que vous vous exposez à un risque de choc électrique, de brûlure, et de blessure avec la lumière du laser.

ATTENTION



Évitez tout risque de choc électrique.

Même si l'instrument a été débranché de toute source d'alimentation électrique, les condensateurs, qui peuvent demeurer chargés jusqu'à 30 secondes, peuvent provoquer un choc électrique.

ATTENTION



Évitez tout risque de brûlure.

La température des composants internes, en particulier celle de la source infrarouge, peut augmenter considérablement pendant le fonctionnement normal de l'instrument. Mettez l'instrument hors tension et attendez 10 minutes au minimum avant de remplacer un de ses composant.

AVERTISSEMENT

Évitez toute lésion corporelle.



- Ne regardez jamais le faisceau laser ni sa réflexion. Ne manipulez ou ne modifiez jamais le laser. Vous pourriez vous exposer à la lumière du laser ou à une haute tension.
- Le réglage du laser, ou toutes procédures non décrites dans les guides et les manuels d'utilisation, pourrait vous exposer à des radiations dangereuses.

REMPLACEMENT DU DESSÉCHANT

Lorsque vous ouvrez le compartiment à desséchant, vous devez empêcher les liquides ou les gaz inflammables d'y pénétrer. Des instructions sur le remplacement du desséchant sont disponibles à la rubrique "Remplacement du desséchant".

DANGER



Évitez tout risque d'explosion.

Avant d'ouvrir le compartiment à desséchant, mettez l'instrument hors tension, débranchez le cordon d'alimentation, et enlevez tous les accessoires et les échantillons du système. L'entrée de liquides ou de gaz inflammables dans le compartiment à desséchant présente un danger d'explosion. Si cela se produit, contactez-nous immédiatement et ne mettez pas l'instrument sous tension tant que le problème n'est pas réglé.

REMPLACEMENT DE LA SOURCE IR

La source IR devient extrêmement chaude lorsque l'instrument fonctionne. Si vous devez remplacer la source IR, évitez les risques de brûlure et d'explosion.

Thermo Scientific

ATTENTION



DANGER

Évitez tout risque d'explosion.

Évitez tout risque de brûlure.



Avant de retirer la source du spectromètre, mettez l'instrument hors tension, débranchez le cordon d'alimentation, débranchez les lignes de purge, et retirez tous les accessoires et échantillons du système. L'entrée de liquides ou de gaz inflammables dans le compartiment de la source peut provoquer une explosion. Si cela se produit, contactez-nous immédiatement et ne mettez pas l'instrument sous tension tant que le problème n'est pas réglé.

La source devient extrêmement chaude pendant le fonctionnement normal du spectromètre. Laissez toujours la source refroidir pendant 10 minutes au minimum

après l'arrêt du spectromètre avant de travailler sur la source.

Aperçu

Le spectromètre Nicolet Summit LITE est un instrument sûr et robuste, mais son utilisation et sa maintenance peuvent comporter des risques. Dans le cadre du fonctionnement normal de l'instrument, faites preuve de prudence lorsque vous manipulez des échantillons et des solvants potentiellement dangereux, et ne touchez pas les parties du spectromètre chaudes ou qui génèrent de la chaleur. Pendant une tâche de maintenance, prenez des précautions pour éviter les blessures initiées par la purge de l'instrument, la manipulation des composants internes ou le remplacement du desséchant.

Déballage et première utilisation du spectromètre

Le déballage et l'installation de votre spectromètre Thermo Scientific Nicolet Summit LITE requiert la mise en place de l'instrument dans votre espace de travail, le raccordement électrique et la mise sous tension de l'instrument, et l'exécution d'un court test de performances.

Avant de commencer

À NOTER

Attendez 24 heures avant de sortir le spectromètre de son sac d'expédition en plastique pour éviter toute formation de condensation à l'intérieur de l'instrument, qui risquerait d'endommager les composants optiques internes. La garantie ne couvre pas les dommages résultant de l'ouverture du sac en plastique scellé avant que l'instrument n'ait atteint la température ambiante.

Rappel : Passez en revue la rubrique Informations relatives à la sécurité et au site d'installation pour vous assurer que votre espace de travail est sûr et adapté à l'instrument.

Déballage et installation

Les étapes de déballage et d'installation de votre spectromètre varient quelque peu selon les accessoires en option que vous avez achetés.

1. Ouvrez la boîte d'expédition

Sortez le spectromètre de la boîte et du sac d'expédition, et retirez tous les matériaux d'emballage supplémentaires.

À NOTER

Soulevez le spectromètre avec précaution.

• Ne touchez pas les fenêtres du compartiment échantillon lors du déballage de l'instrument. Le fait de toucher les fenêtres pourrait provoquer des dommages permanents.



2. (Facultatif) Remplacer les fenêtres du compartiment échantillon.

Si vous avez acheté les fenêtres ZnSe en option, mettez-les en place. Des instructions sur le remplacement des fenêtres sont disponibles à la rubrique <u>"Remplacement des fenêtres du</u> compartiment à échantillons".

- 3. Mise sous tension du spectromètre.
 - a. Raccordez le cordon d'alimentation électrique au spectromètre et à une source de courant alternatif correctement mise à la terre.
 - b. Appuyez sur le bouton d'alimentation de l'instrument pour mettre le spectromètre sous tension (l'allumer). La Barre lumineuse du spectromètre émet une lumière verte cyclique pendant que l'instrument se réchauffe, et affiche une lumière verte continue lorsque l'instrument est prêt à être utilisé.



- 4. Installez le logiciel OMNIC Paradigm.
 - a. Insérez le DVD du logiciel OMNIC Paradigm dans le lecteur.
 - b. Si le DVD ne démarre pas automatiquement, ouvrez les fichiers du DVD et cliquez sur Démarrer pour commencer l'installation. Suivez les instructions à l'écran pour installer le logiciel

OMNIC Paradigm.

L'installation du logiciel peut prendre jusqu'à 20 minutes. Vous devrez redémarrer l'ordinateur pour achever l'installation.

- 5. Connectez le spectromètre
 - a. Utilisez le câble USB Type A/Type B pour raccorder le spectromètre à votre ordinateur. Le spectromètre a une prise USB de Type B.
 - b. Ouvrez le logiciel OMNIC Paradigm. La première fois que vous ouvrez l'application, elle s'ouvre en mode Poste de travail.

🗶 OMNIC	: Paradigm								
		View / Display						💭 Workstation	
Open		le Settings	Deshboard	Spectra Abs /	ST Advanced A	IR Search	Check Quant	y Add To Library	
Measur	ements 🗸 Past w	eek ~			_	Search	0 Ø		Mare
Measuren	sent Name	Las	Modified	lype					
									-
Worldfo	ws ~								More
Name		Date	Created	Last Modified			Preview		
Accessory	Performance Test	8/10	2020 12:00 PM	8/10/2020 12:00					
ATR Acces	nory - PHEUR Qualificatio	n 8/10,	2020 12:00 PM	8/10/2020 12:00	0	- +ł	orn	20	
Nicolet FT	1R - CP Qualification	8/10	2020 12:00 PM	8/10/2020 12:00	1	u		U	
Nicolet FT	1R - Factory Qualification	8/10	2020 12:00 PM	8/10/2020 12:00		9	niant	itic	
Nicolet FT	1R - JP Qualification	8/10	2020 12:00 PM	8/10/2020 12:00	0	0		.mo	
Nicolet FT	IR - PHEUR Qualification	8/10	2020 12:00 PM	8/10/2020 12:00	i .				

c. Une fois que vous avez connecté le spectromètre, le logiciel OMNIC Paradigm le reconnaît automatiquement et affiche le type de l'instrument dans l'angle supérieur droit de l'écran. Si l'instrument ne se connecte pas automatiquement, vous pouvez devoir le faire manuellement.

Pour le connecter manuellement, allez à Configurer > Connectivité et sélectionnez Summit LITE dans la boîte de dialogue de connectivité.

Le logiciel peut employer plusieurs minutes pour démarrer et se connecter à l'instrument. Le logiciel affiche l'état de la connexion dans l'angle supérieur droit de l'écran. Lorsqu'il indique que vous êtes connecté, passez à l'étape suivante.

Note Si vous ouvrez des menus dans le logiciel avant la connexion de l'instrument, les options Aligner et Calibrer seront désactivées. Pour les activer, lorsque l'instrument est connecté, cliquez sur Tableau de bord dans la barre d'outils.

6. Vérification des performances du spectromètre.

Avant de réaliser les étapes suivantes, assurez-vous de l'absence d'accessoires dans le compartiment échantillon.

- a. Alignez les composants optiques du spectromètre.
 - i. Dans le tableau de bord du logiciel OMNIC Paradigm, allez à Acquérir des données
 > Diagnostics > Aligner le spectromètre.
 - ii. Sélectionnez **Démarrer** pour commencer l'alignement. L'écran indique « Alignement terminé » une fois l'alignement terminé avec succès.
- b. Calibrez le laser.
 - i. Sélectionnez Acquérir des données > Diagnostic > Calibration du laser.
 - ii. Sélectionnez Démarrer pour commencer la calibration.

La calibration du laser peut prendre plusieurs minutes. Une fois la calibration terminée, l'écran affiche « Mesure terminée » et indique tout changement de fréquence du laser.

L'instrument peut mettre jusqu'à 12 heures pour atteindre l'équilibre thermique. Pour obtenir les meilleurs résultats, ré-exécutez l'alignement et la calibration après les 12 premières heures de fonctionnement.

- c. Exécutez le workflow Nicolet FTIR Qualification d'usine. Les workflows de vérification des performances et de qualification utilisent du polystyrène comme matériau de référence dans le spectromètre pour exécuter une série de tests standard pour vérifier le fonctionnement de votre instrument et assurer la précision de vos données.
 - i. Revenez au Tableau de bord.
 - ii. Faites défiler le panneau Workflows et sélectionnez Nicolet FTIR Qualification d'usine.
 - iii. Cliquez avec le bouton droit sur le workflow et sélectionnez **Exécuter** pour lancer le workflow.
- 7. Insérez un accessoire d'échantillonnage.

Pour installer l'accessoire de transmission iD1 ou Everest ATR en option, saisissez l'accessoire par les poignées avant et arrière et abaissez l'accessoire pour l'insérer dans le spectromètre. Les accessoires s'insèrent sur deux broches d'alignement de la plaque de base du spectromètre et sont maintenus en place par des aimants puissants.

Première utilisation du spectromètre Nicolet Summit LITE



 Les détails de l'installation d'un accessoire sont disponibles à la rubrique <u>"Installation d'un acces</u>soire d'échantillonnage".

Étapes suivantes

Félicitations ! Votre instrument est prêt à l'emploi. Pour en savoir plus sur l'identification et l'analyse des échantillons avec votre spectromètre, reportez-vous aux guides et tutoriels d'OMNIC Paradigm sur thermofisher.com/ftir-help.

Installation d'un accessoire d'échantillonnage

Le spectromètre Nicolet Summit LITE peut accueillir une grande variété d'accessoires, y compris des accessoires intégrés et d'autres requérant des adaptateurs supplémentaires.

Les accessoires intégrés et les plaques de base sont faciles à installer et à retirer, et sont reconnus automatiquement par le logiciel OMNIC Paradigm.

Installation des accessoires intégrés

Les accessoires intégrés, tels que l'accessoire Everest[™] ATR de Thermo Scientific[™], reposent sur deux broches d'alignement situées sur la plaque de base du spectromètre, et sont maintenus en place par des aimants.



Figure 2-2 : Emplacement des broches d'alignement et des aimants sur un Nicolet Summit.

- 1. Broches
- 2. Aimants
- 3. Trous taraudés

Installer un accessoire intégré

1. Tenez l'accessoire par les poignées avant et arrière et abaissez-le dans le compartiment à échantillons du spectromètre. Des aimants maintiennent l'accessoire en place. Figure 2-3 : Installation d'un accessoire en l'abaissant sur un Nicolet Summit PRO.



2. Si vous purgez votre instrument, raccordez la ligne de purge étiquetée « Vers l'accessoire » à l'entrée de purge à l'arrière de l'accessoire. De plus amples informations sont disponibles à la rubrique <u>"Installation et entretien d'un kit de purge</u>".



Note Pour retirer un accessoire, débranchez la ligne de purge si nécessaire, puis saisissez l'accessoire par les poignées et soulevez-le. Si vous ne comptez pas utiliser un accessoire, stockez-le dans un environnement exempt de poussière, tel qu'une armoire ou une boîte.

Thermo Scientific

Installation des accessoires montés sur la plaque de base

La plaque Adaptateur iD Base permet d'accueillir une grande variété d'accessoires pour l'échantillonnage FTIR, notamment des accessoires de transmission de grande taille et d'autres accessoires spécialisés.



Note Évitez d'installer la plaque Adaptateur iD Base si aucun accessoire n'y est raccordé. Des aimants très puissants maintiennent la plaque en place ; vous devrez peut-être utiliser un outil pour la dégager.

Pour les accessoires montés sur support coulissant, de taille trop importante pour l'accessoire iD1 Transmission, un support coulissant qui s'adapte à l'iD Base est disponible.



La plaque Adaptateur iD Base comporte deux groupes de trous de vis, afin que le support coulissant puisse être installé dans les deux sens. Pour installer le support coulissant, placez-le au-dessus des trous de vis, puis insérez et serrez les deux rondelles et les vis.

Utilisez la vis de réglage de la hauteur pour soulever ou abaisser un échantillon ou un accessoire afin de l'aligner avec le faisceau infrarouge.



Thermo Scientific

- Installer un accessoire à l'aide de la plaque Adaptateur iD Base
- 1. Fixez l'accessoire à la plaque Adaptateur iD Base.
- 2. Placez la plaque Adaptateur iD Base sur les deux broches de la plaque de base du spectromètre. Des aimants puissants maintiennent la plaque et l'accessoire en place.
- 3. Certains accessoires spécialisés montés sur la plaque de base sont fixés à l'instrument en utilisant 4 vis fournies avec l'accessoire.

Installation des accessoires montés sur support coulissant

L'accessoire iD1 Transmission est conçu pour accueillir des cellules de transmission de gaz ou de liquide et des supports de films minces ou de pastille qui ont un support coulissant de 2 x 3 po.



- Installer un accessoire monté sur support coulissant
- 1. Faites glisser l'accessoire dans deux fentes de l'accessoire iD1 Transmission.

Choisissez les deux fentes qui aligneront le centre de l'accessoire avec le point focal du faisceau



(désigné par des flèches sur le boîtier de l'accessoire).

Thermo Scientific

Comprendre la barre lumineuse du spectromètre

Le spectromètre Thermo Scientific Nicolet Summit LITE est équipé d'une barre lumineuse pratique, destinée à fournir une indication visuelle rapide (en un coup d'œil) de la qualité du produit et de l'état de l'instrument.

Les tableaux suivants décrivent les significations des différents signaux de la barre lumineuse.

Signal	État du système	Description
Lumière verte cyclique	Le système est sous tension et en cours de stabilisation thermique.	La lumière verte se déplace de gauche à droite de manière répétée. La durée de stabilisation thermique est d'environ 2 minutes.

STABILISATION THERMIQUE EN COURS.

ACQUISITION DE DONNÉES TRADITIONNELLE

Signal	État du système	Description
Lumière verte continue	Prêt à l'emploi	La barre lumineuse entière émet une lumière verte continue

Première utilisation du spectromètre Nicolet Summit LITE



ACQUISITION DE DONNÉES AVEC LA FONCTIONNALITÉ SMART BACKGROUND

Signal	État du système	Description
Pulsation bleue	Acquisition d'un Smart Background	La barre lumineuse entière émet des pulsations bleues
Lumière bleue passante	Acquisition de données d'échantillon (initiée par l'utilisateur)	Une petite section de lumière bleue effectue des allers-retours

RÉSULTATS DE L'ANALYSE

Signal	État du système	Description
Lumière verte, avec un % de remplissage	Pourcentage d'identification ou résultat du QCheck supérieur au seuil	Remplissage en vert au prorata du pourcentage d'identification. Par exemple, environ 90 % de remplissage pour un pourcentage d'identification de 90.
Lumière orange avec un % de remplissage	Pourcentage d'identification ou résultat du QCheck inférieur au seuil	Remplissage en orange au prorata du pourcentage d'identification. Par exemple, environ 30 % de remplissage pour un pourcentage d'identification de 30.
DIAGNOSTICS		
Signal	État du système	Description
Lumière rouge clignotante	Erreur système	La barre lumineuse entière clignote en rouge. Une description de l'erreur est disponible à la rubrique État du système du logiciel OMNIC Paradigm.

Consultation de vos données via l'application OMNIC Anywhere

OMNIC Anywhere est une application reposant sur le cloud qui vous permet de visualiser, d'examiner ou de partager vos données depuis un PC, un ordinateur Apple, un périphérique Android ou iOS connecté.

Figure 2-4 : OMNIC Anywhere dans le navigateur web



Un spectromètre Nicolet Summit LITE conjointement à un compte Connect gratuit, vous permet de mesurer des échantillons depuis une salle de classe ou un laboratoire, de télécharger les données sur votre compte Connect, et de visualiser, d'examiner ou de partager les données avec un autre appareil situé dans un espace privé ou un espace de travail.

Des instructions sur la création d'un compte Connect et la consultation de vos données à l'aide de l'application OMNIC Anywhere sont disponibles à la rubrique « <u>Spectroscopie FTIR activée sur le</u> <u>cloud</u> ».

Chapitre 3

Options et accessoires

Le spectromètre Nicolet Summit LITE peut être utilisé avec le kit de purge en option, et est compatible avec une large gamme d'accessoires d'échantillonnage.

Accessoires d'échantillonnage pour le spectromètre Summit	32
Installation et entretien d'un kit de purge	35

Accessoires d'échantillonnage pour le spectromètre Summit

Les spectromètres Thermo Scientific Nicolet Summit LITE FTIR sont compatibles avec les accessoires d'échantillonnage Thermo Scientific[™] Nicolet[™] iS5 iD et des centaines d'autres accessoires de tiers.

Des instructions sur l'installation d'un accessoire d'échantillonnage sont disponibles à la rubrique "Accessoires d'échantillonnage pour le spectromètre Summit".

Grâce aux accessoires et adaptateurs suivants, le spectromètre Summit peut mesurer rapidement et facilement une vaste gamme d'échantillons.

ACCESSOIRE EVEREST ATR



L'accessoire Everest[™] ATR est un accessoire de réflectance totale atténuée à une seule réflexion, polyvalent et performant, idéal pour l'analyse de liquides, de solides, de pâtes ou de poudres. L'accessoire Everest ATR a été spécialement conçu pour s'adapter aux spectromètres Summit.

Contactez-nous pour en savoir plus.

ACCESSOIRE ID1 TRANSMISSION



L'accessoire iD1 Transmission est pratique pour l'échantillonnage de substances sur films, en granulés, de liquides et gazeuses, et s'adapte à une vaste gamme d'accessoires de transmission. Le compartiment à échantillons peut accueillir des cuvettes et des cellules à gaz mesurant jusqu'à 10 cm.

Contactez-nous pour en savoir plus.

ACCESSOIRES ID5/ID7 ATR



Les accessoires iD5 et iD7 ATR (conçus pour le spectromètre Nicolet[™] iS5) peuvent être utilisés comme une alternative à l'accessoire Everest ATR. Cependant, bien que les accessoires ATR soient d'excellents accessoires polyvalents, l'utilisation de l'accessoire Everest ATR avec les spectromètres Summit reste recommandée.

De plus amples informations sont disponibles aux rubriques <u>« Accessoire iD5 ATR pour spectromètre</u> Nicolet[™] iS5 » ou « Accessoire iD7 ATR pour spectromètre Nicolet[™] iS5 ».

PLAQUE ADAPTATEUR ID BASE



La plaque Adaptateur iD Base permet d'accueillir une grande variété d'accessoires tiers pour l'échantillonnage FTIR, y compris des accessoires de transmission de grande taille et d'autres accessoires spécialisés. Associez la plaque Adaptateur iD Base avec le porte-échantillons à support coulissant pour obtenir une plus grande flexibilité.

De plus amples informations sont disponibles à la rubrique « Adaptateur iD Base ».

PORTE-ÉCHANTILLON À SUPPORT COULISSANT



Associez un porte-échantillon à support coulissant avec la plaque Adaptateur iD Base pour pouvoir

utiliser les accessoires à support coulissant qui sont trop grands pour l'accessoire iD1 Transmission.

Contactez-nous pour en savoir plus.

ACCESSOIRE GOLDEN GATE ATR



L'accessoire Golden Gate ATR, qui comporte une "enclume" en saphir pour les contacts haute pression et est doté d'un boîtier robuste, vous permet d'analyser toute une gamme d'échantillons - des fibres et particules simples ou des solides durs aux liquides corrosifs. L'accessoire Golden Gate ATR est également idéal pour le macro-échantillonnage.

De plus amples informations sont disponibles à la rubrique <u>« Thermo Scientific™ Golden Gate ATR pour</u> spectromètre Nicolet iS5 FTIR ».

Installation et entretien d'un kit de purge

La purge du spectromètre avec de l'air sec ou de l'azote protège les composants internes de l'humidité et d'autres contaminants environnementaux. Pour installer le kit de purge, vous devez assembler les vannes et les régulateurs, remplacer le support de desséchant, raccorder le gaz de purge et régler sa pression et son débit.

AVERTISSEMENT



Évitez tout risque d'explosion.

N'utilisez jamais de gaz inflammable, combustible ou toxique pour purger cet instrument. Le gaz de purge ne doit pas contenir d'huile ni d'autres matières réactives. La chaleur provenant de la source ou de l'absorption du laser peut enflammer des gaz inflammables ou des matières réactives dans le gaz de purge. Utilisez uniquement de l'azote ou de l'air pur et sec pour purger l'instrument.

À NOTER

Nous recommandons de garder le spectromètre scellé et rempli de desséchant et/ou purgé à tout moment. Les dommages causés à l'instrument par un défaut d'étanchéité et de desséchant et/ou de purge ne sont pas couverts par la garantie. Si vous avez des questions concernant cette exigence, contactez-nous.

CONFIGURATION REQUISE

Avant d'installer le kit de purge, vous aurez besoin d'une source d'air sec ou d'azote répondant aux spécifications Thermo Scientific pour le spectromètre Nicolet™ Summit. De plus amples informations sont disponibles dans le Guide de sécurité et d'installation sur site de l'instrument.

Note Pour de meilleurs résultats, asséchez le gaz de purge jusqu'à atteindre un point de rosée inférieur ou égal à -70 °C (-94 °F).

MATÉRIEL REQUIS

Votre kit de purge comprend les éléments suivants :

- Collecteur de purge à deux zones
- Support de desséchant de purge, avec cartouche de desséchant de purge et joint torique

En plus du kit de purge, vous aurez besoin des outils et du matériel suivants :

- Une clé anglaise de 3/4 po
- Une clé anglaise de 11/16 po
- Un tournevis cruciforme n° 2
- Du ruban d'étanchéité pour filetage (« ruban de plombier » ou « ruban téflon »)

Installer un kit de purge

- 1. Fixez l'ensemble de tuyauterie à la source du gaz de purge.
 - a. Installez la vanne de régulation et un raccord mâle de 1/4" ou un raccord femelle de 3/8" sur la source du gaz de purge (choisissez une vanne et des raccords qui conviennent à la source du gaz de purge).



b. Si vous avez installé un raccord mâle de 1/4 po, passez à l'étape suivante.

Si vous avez installé un raccord femelle de 3/8" sur la source du gaz de purge, installez le manchon de réduction de 3/8" à 1/4" inclus dans votre kit de purge. Enveloppez le manchon de réduction avec du ruban d'étanchéité pour filetage avant de l'installer, et serrez le raccord à l'aide d'une clé de 11/16".



c. Enveloppez le manchon de réduction ou le raccord mâle de 1/4 po avec du ruban d'étanchéité pour filetage, puis installez le raccord de pression. Serrez le raccord à l'aide d'une clé anglaise de 3/4".



d. Emboîtez fermement l'entrée mâle de l'ensemble de tuyauterie murale dans le raccord à dégagement rapide.



- 1. Vers l'accessoire
- 2. Vers le spectromètre
- 3. Débitmètre
- 4. Régulateur de pression
- 5. Raccord à déclenchement rapide
- 6. Ensemble de tuyauterie murale
- 2. Remplacez la support de desséchant.
 - a. Mettez l'instrument hors tension, débranchez le cordon d'alimentation et retirez tous les accessoires et échantillons du système.

AVERTISSEMENT

Évitez tout risque d'explosion.



Avant d'ouvrir le compartiment à desséchant, mettez l'instrument hors tension, débranchez le cordon d'alimentation et retirez tous les accessoires et les échantillons du système. L'entrée de liquides ou de gaz inflammables dans le compartiment à desséchant présente un danger d'explosion. Si cela se produit, contactez-nous immédiatement et ne mettez pas l'instrument sous tension tant que le problème n'est pas réglé.

Thermo Scientific

b. Desserrez les deux vis à tête cruciforme captives de support de desséchant standard, et soulevez le support verticalement pour le sortir de l'instrument.



Note Si le support de desséchant standard est stocké dans une boîte sèche ou un sac scellé (vous pouvez utiliser le sac dans lequel le support de purge a été livré), le support peut être réutilisé.

c. Retirez et jetez le grand joint torique.



- d. Ouvrez le kit de purge et le sachet de desséchant scellé.
- e. Retirez le joint torique du sachet et installez-le sur la plaque de base de l'instrument (voir l'image précédente). Appuyez sur le nouveau joint torique (au lieu de le rouler) pour le mettre en place, en veillant à ce qu'il soit bien logé dans l'encoche.
- f. Retirez le support de desséchant avec les raccords de purge du paquet, insérez le support dans le compartiment à desséchant, assurez-vous qu'il est correctement positionné sur le joint torique, puis fixez-le à l'aide des deux vis.

Vérifiez que l'orientation du support est correct. Lorsque le support est installé, vous devez être en mesure de lire l'inscription sur l'indicateur d'humidité depuis l'avant de l'instrument.



- 1. Entrée de purge
- 2. Vanne de décharge
- 3. Purgez le support de desséchant.
- 3. Connectez la source du gaz de purge à l'instrument.
 - a. Raccordez le coupleur de débit de l'ensemble de tuyauterie murale (étiqueté « Vers le spectromètre ») à l'entrée de purge du support de desséchant.
 - b. Remettez en place tout accessoire d'échantillonnage que vous avez retiré précédemment.
 - c. Si vous purgez l'accessoire, raccordez le coupleur de débit droit de l'ensemble de tuyauterie (étiqueté « Vers l'accessoire ») à l'entrée de purge de l'accessoire installé.
- 4. Réglez les commandes de gaz de purge.
 - a. Ouvrez la vanne de régulation principale et réglez les commandes de gaz de purge comme indiqué ci-dessous (de plus amples informations sont disponibles à la rubrique Réglage des commandes de gaz de purge). Si vous ne purgez pas l'accessoire, réglez le régulateur de pression des accessoires sur zéro.

Matériel	Pression (psig)	Plage de débit (scfh)
Spectromètre Summit	5	1

Matériel	Pression (psig)	Plage de débit (scfh)
Accessoire Everest	5	4-10
Accessoire iD1 Transmission	5	4-10
Accessoire iD3 ATR	5	4-10
Accessoire iD5 ATR	5	4-10
Accessoire iD7 ATR	5	4-10
iD Foundation	5	4-10

- b. Raccordez le cordon d'alimentation à l'instrument et mettez ce dernier sous tension.
- c. Attendez 30 à 60 minutes pour que l'instrument soit complètement purgé.

Réglage des commandes de gaz de purge

En réglant correctement les commandes de gaz de purge, vous protégerez votre spectromètre de l'humidité, sans générer de vibrations.

Pour de meilleurs résultats, asséchez le gaz de purge jusqu'à atteindre un point de rosée de -70 °C (-94 °F) ou inférieur.

- Régler les commandes de gaz de purge
- 1. Ouvrez la vanne principale pour démarrer l'écoulement du gaz de purge dans le régulateur.
- 2. Ajustez le régulateur de pression du spectromètre jusqu'à ce que le manomètre indique une pression de 5 psig (34 kPa).



- 1. Réglage du débit de l'accessoire
- 2. Réglage de la pression de l'accessoire
- 3. Réglage du débit du spectromètre
- 4. Réglage de la pression du spectromètre
- 3. Réglez le débitmètre du spectromètre sur 1 scfh (0,47 l/min).
- 4. Si vous utilisez un accessoire qui est purgé, réglez les commandes du gaz de purge pour l'accessoire comme indiqué ci-dessous.
 - Pression de l'accessoire : 5 psig (34 kPa)
 - Débit de l'accessoire : 4 à 10 scfh (1,9 à 4,7 l/min)

À NOTER

Des débits supérieurs aux valeurs recommandées peuvent provoquer des vibrations susceptibles d'affecter la qualité de vos données.

Vérification et remplacement du filtre de gaz de purge

Remplacez le filtre lorsqu'il devient jaune, ou présente une quelconque décoloration, ou s'il est contaminé par des débris ou des particules étrangères.

À NOTER

Nous recommandons de garder constamment le spectromètre scellé, dessiqué et/ou purgé. Les dommages causés à l'équipement par un défaut d'étanchéité et de desséchant et/ou de purge ne sont pas couverts par la garantie. Si vous avez des questions concernant cette exigence, contactez-nous.

Figure 3-1 : Le filtre de purge est situé à l'intérieur de la cuve en plastique, sous le manomètre



Note Pour toute commande de pièces, contactez le support technique.

Remplacer le filtre de gaz de purge

- 1. Fermez le débit du gaz de purge au niveau de la vanne principale. Ne baissez pas le débitmètre ni le régulateur de pression.
- 2. Retirez la cuve en plastique qui abrite le filtre, puis retirez le filtre (vous pouvez les dévisser tous les deux à la main).



3. Installez le nouveau filtre, puis réinstallez la cuve.

4. Ouvrez la vanne principale pour activer le débit de purge vers l'instrument, et vérifiez le débit correct du gaz de purge.

Chapitre 4

Maintenance

Votre spectromètre Nicolet Summit LITE FTIR nécessite peu d'entretien, et aucune maintenance quotidienne. Cependant, quelques soins de base permettront d'assurer les performances optimales régulières de l'instrument.

Calendrier de maintenance du spectromètre Summit	46
Alignement des composants optiques du spectromètre	50
Calibration de la fréquence du laser	51
Nettoyage du spectromètre	52
Remplacement du desséchant	53
Remplacement de la source	59
Remplacement des fenêtres du compartiment à échantillons	63

Calendrier de maintenance du spectromètre Summit

Votre spectromètre Thermo Scientific Nicolet Summit LITE FTIR nécessite peu d'entretien, et aucune maintenance quotidienne. Cependant, quelques soins de base permettront d'assurer les performances optimales régulières de l'instrument. Suivez les directives de cette rubrique. Pour obtenir les meilleures performances, laissez le spectromètre sous tension.

À NOTER

L'électricité statique peut endommager de façon permanente les composants critiques de votre spectromètre. Pour éviter de tels dommages, prenez en compte les recommandations suivantes :

- Avant de débrancher le cordon d'alimentation électrique, déchargez l'électricité statique que vous avez pu accumuler en touchant la base métallique du spectromètre.
- Laissez les pièces de rechange dans leur emballage de protection jusqu'à ce que vous soyez prêt (e) à les installer dans l'instrument.

Entretien hebdomadaire

VÉRIFICATION DES PERFORMANCES DU SPECTROMÈTRE

Le logiciel OMNIC Paradigmcomprend des workflows de qualification et de vérification des performances (PV) conçus pour votre spectromètre Summit. Les workflows de qualification sont des tests de qualification standard à l'échelle de l'industrie que vous reconnaitrez probablement (par ex., Pharmacopée européenne, Pharmacopée japonaise, etc.). Le workflow PV exécute une série de tests standard destinée à vérifier le fonctionnement de l'instrument et à garantir l'exactitude de vos données. Tous les étalons de référence requis sont fournis à l'intérieur de l'instrument et sont contrôlés par le logiciel OMNIC Paradigm.

Nous vous recommandons d'exécuter le workflow PV ou de qualification de votre choix une fois par semaine au minimum.

NETTOYAGE DU SPECTROMÈTRE

Le spectromètre et l'écran tactile doivent être nettoyés uniquement en suivant les recommandations. Reportez-vous à <u>"Nettoyage du spectromètre"</u>.

Entretien mensuel

VÉRIFICATION DE L'INDICATEUR D'HUMIDITÉ

Les composants optiques du spectromètre peuvent être facilement endommagés par un taux d'humidité excessif dans l'air. Le spectromètre est scellé, et ses composants sont protégés par deux cartouches de desséchant qui absorbent l'humidité. Surveillez l'humidité de l'instrument dans le logiciel OMNIC Paradigm.

Les composants optiques du spectromètre peuvent être facilement endommagés par un taux d'humidité excessif dans l'air. Le spectromètre est scellé, et ses composants sont protégés par deux cartouches de desséchant qui absorbent l'humidité. Un indicateur d'humidité (voir son emplacement dans la figure cidessous) surveille le taux d'humidité à l'intérieur du spectromètre.

Vous pouvez également surveiller l'humidité avce le logiciel OMNIC Paradigm. Dans l'interface Bureau, cliquez sur l'icône de l'instrument pour afficher l'état du système ou sur l'interface de l'écran tactile, sélectionnez l'icône Diagnostics. Le niveau d'humidité est indiqué dans la section Environnement.





Vérifiez l'indicateur d'humidité au minimum une fois par mois, et changez les cartouches de desséchant selon les besoins (voir le tableau ci-dessous), ou achetez et installez un kit de purge. De plus amples informations sont disponibles à la rubrique <u>"Installation et entretien d'un kit de purge"</u>.

État de l'indicateur d'humidité	Signification	Action
Bleu	Le desséchant est conforme	None (Aucune)
Bleu clair	Le desséchant devient saturé d'humidité et n'offre plus une protection suffisante	Remplacez le desséchant.
Rose ou blanc	Le desséchant est expiré.	Remplacez le desséchant et l'indicateur d'humidité

Tableau 1 - États des indicateurs d'humidité et recommandations

De plus amples informations sont disponibles à la rubrique "Remplacement du desséchant".

À NOTER

Nous vous recommandons de constamment garder le spectromètre scellé et dessiqué, ou scellé et purgé avec de l'air sec ou de l'azote. Les dommages causés à l'équipement par un défaut d'étanchéité et de desséchant et/ou de purge ne sont pas couverts par la garantie. Si vous avez des questions concernant cette exigence, contactez-nous.

VÉRIFICATION DU FILTRE DE GAZ DE PURGE

Si votre spectromètre est purgé à l'azote ou à l'air sec, vérifiez le filtre de purge une fois par mois au minimum.

Remplacez le filtre lorsqu'il devient jaune, ou présente une quelconque décoloration, ou s'il est contaminé par des débris ou des particules étrangères. De plus amples informations sont disponibles à la rubrique <u>"Vérification et remplacement du filtre de gaz de purge"</u>.

Figure 4-2 : Le filtre de purge est situé à l'intérieur de la cuve en plastique, sous le manomètre



Alignement des composants optiques du spectromètre

Alignez les composants optiques de votre spectromètre si un worflow de qualification, tel que le test Nicolet FTIR - Qualification d'usine, indique un échec. L'alignement du spectromètre optimise l'énergie qui atteint le détecteur et maximise le signal du détecteur. Si un test de performance ou de qualification échoue, alignez les composants optiques du spectromètre, calibrez la fréquence du laser et refaites le test. Si le test échoue à nouveau, contactez un représentant du service technique local pour obtenir de l'aide.

- Aligner les composants optiques du spectromètre (interface écran tactile)
- 1. Dans l'écran d'accueil, sélectionnez l'icône Diagnostics [
- 2. Ouvrez l'onglet Aligner et sélectionnez Aligner.

Le système aligne automatiquement le spectromètre et indique la fin de l'alignement.

- Aligner les composants optiques du spectromètre (interface opérateur)
- 1. Naviguez jusqu'à la zone Administrateur et sélectionnez l'icône Diagnostics [
- 2. Ouvrez l'onglet Aligner et sélectionnez Aligner.

Le système aligne automatiquement le spectromètre et indique la fin de l'alignement.

- Aligner les composants optiques du spectromètre (interface bureau)
- 1. Sélectionnez Acquérir des données > Diagnostics > Aligner le spectromètre.
- 2. Dans la boîte de dialogue Aligner le spectromètre, cliquez sur Démarrer.

Le système aligne automatiquement le spectromètre et indique la fin de l'alignement.

Calibration de la fréquence du laser

Calibtez la fréquence du laser si un workflow de qualification, tel que le workflow Nicolet FTIR -Qualification d'usine, indique que le test a échoué. Lorsqu'un test échoue, il convient d'abord d'aligner le spectromètre, puis de calibrer la fréquence du laser et de recommencer le test. Si le test échoue à nouveau, contactez un représentant du service technique local pour obtenir de l'aide.

Calibrer le laser (interface écran tactile)

- 1. Dans l'écran d'accueil, sélectionnez l'icône Diagnostics [
- 2. Ouvrez l'onglet Calibration du laser et sélectionnez Calibrer pour commencer l'opération.

La calibration du laser terminée, un message indique si la fréquence du laser a été modifiée, et affiche la nouvelle fréquence.

Calibrer le laser (interface opérateur)

- 1. Naviguez jusqu'à la zone Administrateur et sélectionnez l'icône Diagnostics [1] pour ouvrir la vue Diagnostics.
- 2. Ouvrez l'onglet Calibration du laser et sélectionnez Calibrer pour commencer l'opération.

La calibration du laser terminée, un message indique si la fréquence du laser a été modifiée, et affiche la nouvelle fréquence.

Calibrer le laser (interface bureau)

- 1. Sélectionnez Acquérir des données > Diagnostic > Calibration du laser.
- 2. Dans la boîte de dialogue Calibration du laser, cliquez sur Démarrer pour calibrer le laser.

Lorsque le système a terminé la calibration du laser, un message indique si la fréquence a été modifiée, et affiche la nouvelle fréquence.

Nettoyage du spectromètre

Avant de nettoyer le spectromètre Thermo Scientific Nicolet Summit LITE, mettez-le hors tension et débranchez le cordon d'alimentation.

Nettoyage du spectromètre Summit

Nettoyez délicatement l'extérieur de l'instrument avec un chiffon doux et propre légèrement humidifié avec du savon doux.

De la poussière peut s'accumuler sur le boîtier électronique à l'arrière de l'instrument, au risque d'interférer avec la dissipation de la chaleur, ce qui peut réduire la durée de vie des composants électroniques.

Pour enlever la poussière accumulée à l'arrière de l'instrument, utilisez de l'air comprimé pour la chasser. N'utilisez pas de liquide pour enlever la poussière à l'arrière du spectromètre.

ATTENTION



Évitez tout risque de choc électrique. Pour éviter tout risque de choc électrique, ne laissez jamais du liquide s'écouler dans l'alimentation électrique ou dans la partie arrière de l'instrument.

À NOTER

N'utilisez pas de détergents agressifs, de solvants, de produits chimiques ou d'abrasifs, sous peine d'endommager la finition du système. Ne laissez pas un liquide entrer en contact avec les fenêtres d'isolation du compartiment à échantillons.

À NOTER

Les fenêtres peuvent être rayées et abîmées très facilement. Ne les touchez pas et n'essayez pas de les nettoyer. La poussière n'affectera pas le signal, tandis que les traces de doigt peuvent dégrader les performances de l'instrument et endommager de façon permanente les miroirs ou les fenêtres. Si vous souhaitez enlever la poussière à la surface d'un miroir ou d'une fenêtre, chassez-la avec un jet d'air ou d'azote propre et sec à basse pression. Ne laissez jamais un liquide entrer en contact avec une fenêtre d'isolation ou un composant optique de l'instrument.

Remplacement du desséchant

Si votre instrument n'est pas équipé d'un kit de purge, ou si la purge est désactivée, vous devez surveiller le niveau d'humidité à l'intérieur du spectromètre. Remplacez le desséchant lorsque le papier indicateur d'humidité devient rose (il peut être rose clair ou presque blanc). Remplacez le papier indicateur d'humidité chaque fois que vous remplacez le desséchant.

Pour remplacer le desséchant, vous devez retirer les cartouches de desséchant et les remplacer ainsi que le joint torique.

OUTILS REQUIS

Avant de commencer, assurez-vous de réunir les éléments suivants :

- Un kit de remplacement du desséchant
- Une clé hexagonale de 0,05 po
- Un tournevis cruciforme n° 2
- Des gants, doigtiers ou lingettes de laboratoire (pour manipuler l'indicateur d'humidité).

AVERTISSEMENT

Évitez tout risque d'explosion.



Avant d'ouvrir le compartiment à desséchant, mettez l'instrument hors tension, débranchez le cordon d'alimentation et retirez tous les accessoires et les échantillons du système. L'entrée de liquides ou de gaz inflammables dans le compartiment à desséchant présente un danger d'explosion. Si cela se produit, contactez-nous immédiatement et ne mettez pas l'instrument sous tension tant que le problème n'est pas réglé.

À NOTER

Assurez-vous que rien ne tombe dans l'instrument lorsque le couvercle du desséchant est retiré.

AVERTISSEMENT



Remplacez le desséchant uniquement avec les pièces de rechange que nous fournissons.

Remplacer le desséchant

- 1. Mise hors tension du spectromètre.
 - a. Pour éteindre le spectromètre, appuyez sur le bouton d'alimentation.
 - b. Débranchez le cordon d'alimentation.
 - c. Arrêter la purge, si applicable, et débranchez la ligne de purge de l'instrument et de tout accessoire.

Note Vous pouvez utiliser un support de desséchant avec ou sans connecteur de purge pour maintenir l'humidité requise. Les illustrations ci-dessous présente une support de desséchant « sans purge ». Les instructions pour le remplacement des cartouches de desséchant sont identiques pour les deux types de support.

- 2. Retrait de la cartouche de desséchant.
 - a. Retirez tout accessoire installé et tous les échantillons de l'instrument.
 - b. Utilisez le tournevis cruciforme n° 2 pour desserrer les deux vis captives du support de desséchant, et soulevez le support verticalement pour l'extraire de l'instrument.



Thermo Scientific

c. Retirez et jetez le grand joint torique.



Note Veillez à ne pas faire tomber le joint torique dans le compartiment à desséchant.

3. Retrait des cartouches de desséchant et remplacement de l'indicateur d'humidité.

AVERTISSEMENT Image: Avertise of the second secon

a. Utilisez la clé hexagonale de 0,05 po pour desserrer (d'environ deux tours) les huit vis qui maintiennent les deux cartouches de desséchant saturé dans le support de desséchant. Retirez les cartouches.

A	X			8			0	7	1
			100					2	
2	Server		01				0 0' 0 0		
	10.0000		G			4 A			
	9		E O	6.6.8 17.4.8	0.0		9 0 9 9	3	
	3		NIO						
	3.1		Ŧ						
	B.s.	••	12	• • •					
		800	-						
	2000000	222	ž	* * *	00	a a	* *	21	
	12	000	-	5.0.0		a a			
	11	000			* *	5 0 5 5	* *		
	S	•• 3	E		• •		-	3	

- 1. Support de desséchant
- 2. Retirez ces vis (quatre de chaque côté)
- 3. Cartouches de desséchant
- b. Retournez le support de desséchant et détachez l'indicateur d'humidité usagé de la fenêtre. Jetez l'indicateur usagé.

À NOTER

Portez toujours des gants de laboratoire ou des doigtiers, ou utilisez une lingette de laboratoire, lorsque vous manipulez un nouvel indicateur d'humidité. La graisse ou l'humidité de la peau peut décolorer l'indicateur.

- 4. Installation des nouvelles cartouches de desséchant.
 - a. Ouvrez le paquet scellé qui contient le nouveau desséchant (qui doit être scellé pour préserver l'intégrité du produit) et retirez le nouvel indicateur d'humidité.
 - b. Retournez le support de desséchant et appuyez l'indicateur bleu sur la fenêtre de sorte que les bords plats de l'indicateur soient alignés avec les bords du support. Le texte doit être visible à

travers la fenêtre lorsque vous retournez le support.

À NOTER

Vérifiez que l'indicateur est bien appuyé contre la fenêtre afin qu'il ne se détache pas dans le compartiment à desséchant lorsque vous réinstallerez le support.

- c. Insérez les cartouches dans le support de desséchant et serrez les huit vis (d'environ 2 tours) jusqu'à ce qu'elles soient au fond du support et à la même hauteur.
- 5. Insertion du support de desséchant.
 - a. Retirez le nouveau joint torique du paquet et installez-le sur la plaque de base de l'instrument.

Appuyez sur le nouveau joint torique (au lieu de le rouler) pour le mettre en place, en veillant à ce qu'il soit bien logé dans la rainure.

b. Insérez le support de desséchant dans le spectromètre afin que le texte de l'indicateur d'humidité puisse être lu par une personne se tenant devant l'instrument.

Assurez-vous que le support est correctement placée sur le joint torique, puis serrer les deux vis captives à l'aide du tournevis cruciforme n° 2.

- 6. Mise sous tension du spectromètre.
 - a. Raccordez le cordon d'alimentation à l'instrument et mettez ce dernier sous tension.
 - b. Rebranchez la ligne de purge sur l'instrument, si applicable, et activez la purge.
 - c. Reconnectez le spectromètre à l'ordinateur.
- 7. Vérification des performances du spectromètre.

Avant de réaliser les étapes suivantes, assurez-vous de l'absence d'accessoires dans le compartiment échantillon.

- a. Alignez les composants optiques du spectromètre.
 - i. Dans le tableau de bord du logiciel OMNIC Paradigm, allez à Acquérir des données
 > Diagnostics > Aligner le spectromètre.
 - ii. Sélectionnez **Démarrer** pour commencer l'alignement. L'écran indique « Alignement terminé » une fois l'alignement terminé avec succès.

- b. Calibrez le laser.
 - i. Sélectionnez Acquérir des données > Diagnostic > Calibration du laser.
 - ii. Sélectionnez Démarrer pour commencer la calibration.

La calibration du laser peut prendre plusieurs minutes. Une fois la calibration terminée, l'écran affiche « Mesure terminée » et indique tout changement de fréquence du laser.

L'instrument peut mettre jusqu'à 12 heures pour atteindre l'équilibre thermique. Pour obtenir les meilleurs résultats, ré-exécutez l'alignement et la calibration après les 12 premières heures de fonctionnement.

- c. Exécutez le workflow Nicolet FTIR Qualification d'usine. Les workflows de vérification des performances et de qualification utilisent du polystyrène comme matériau de référence dans le spectromètre pour exécuter une série de tests standard pour vérifier le fonctionnement de votre instrument et assurer la précision de vos données.
 - i. Revenez au Tableau de bord.
 - ii. Faites défiler le panneau Workflows et sélectionnez Nicolet FTIR Qualification d'usine.
 - Cliquez avec le bouton droit sur le workflow et sélectionnez Exécuter pour lancer le workflow.

Remplacement de la source

La source IR est facilement accessible depuis la partie inférieure du spectromètre, et peut être remplacée sans avoir à retirer le couvercle du spectromètre.

OUTILS ET MATÉRIEL REQUIS

- Un tournevis cruciforme n° 1
- Un kit de remplacement de source

ATTENTION



Avant de remplacer la source, mettez l'instrument hors tension. Remplacez la source uniquement avec les pièces de rechange que nous fournissons.

Remplacer la source

- 1. Mise hors tension et débranchement du spectromètre.
 - a. Mettez l'instrument hors tension et débranchez le cordon d'alimentation.
 - b. Débranchez tous les câbles de l'instrument, tels que le câble Ethernet ou les périphériques USB.
 - c. Si le spectromètre ou un accessoire installé est purgé, débranchez les lignes de purge du spectromètre et de l'accessoire (les raccords rapides arrêtent automatiquement le flux). De plus amples informations sont disponibles à la rubrique "Installation et entretien d'un kit de purge".
 - d. Retirez tous les accessoires installés sur l'instrument.

ATTENTION



Évitez tout risque de brûlure.

La source devient extrêmement chaude lorsque l'instrument est en marche. Attendez 10 minutes au minimum après avoir mis l'instrument hors tension avant de passer à l'étape suivante.

- 2. Retrait de la source.
 - a. Inclinez doucement l'instrument vers l'arrière jusqu'à ce qu'il repose sur le boîtier électronique.



b. Utilisez un tournevis cruciforme n° 1 pour desserrer complètement les trois vis captives qui maintiennent la source en place.



c. Sortez la poignée en la faisant pivoter. Tenez l'instrument d'une main et tirez fermement sur la poignée pour faire sortir la source directement de l'instrument.



À NOTER

Ne touchez jamais un élément de la source avec vos doigts nus. La graisse de la peau ou d'autres dépôts réduiraient sa durée de vie. Utilisez toujours des doigtiers, des gants ou des lingettes de laboratoire propres lorsque vous manipulez un élément de la source.

- 3. Insérez la nouvelle source.
 - a. Assurez-vous que les trois vis captives sont bien droites, puis insérez lentement la source dans la cavité jusqu'à ce que le support de la source repose contre l'instrument.
 - b. Serrez les vis pour fixer la source et mettez la poignée en position de stockage.
- 4. Rebranchez le spectromètre et mettez-le sous tension.
 - a. Remettez soigneusement l'instrument en position verticale et rebranchez les câbles que vous aviez débranchés.
 - b. Remplacez le desséchant. De plus amples informations sont disponibles à la rubrique "Remplacement du desséchant".

Les composants optiques internes étant exposés à l'atmosphère lors du retrait de la source, le desséchant doit toujours être remplacé lorsque la source vient d'être remplacée.

AVERTISSEMENT



Évitez tout risque de choc électrique.

Pour assurer une mise à la terre satisfaisante et éviter les risques de choc électrique, n'utilisez pas une prise de courant reliée à la terre d'un conduit. Le fil de terre doit être un fil ne transportant pas de courant et doit être raccordé à la terre sur le tableau de distribution principal.

c. Raccordez le cordon d'alimentation à l'instrument et appuyez sur le bouton d'alimentation pour mettre le spectromètre sous tension.

Note N'installez aucun accessoire d'échantillonnage avant d'avoir aligné le spectromètre et vérifié ses performances

5. Vérification des performances du spectromètre.

Avant de réaliser les étapes suivantes, assurez-vous de l'absence d'accessoires dans le compartiment échantillon.

- a. Alignez les composants optiques du spectromètre.
 - i. Dans le tableau de bord du logiciel OMNIC Paradigm, allez à Acquérir des données
 > Diagnostics > Aligner le spectromètre.
 - ii. Sélectionnez **Démarrer** pour commencer l'alignement. L'écran indique « Alignement terminé » une fois l'alignement terminé avec succès.
- b. Calibrez le laser.
 - i. Sélectionnez Acquérir des données > Diagnostic > Calibration du laser.
 - ii. Sélectionnez Démarrer pour commencer la calibration.

La calibration du laser peut prendre plusieurs minutes. Une fois la calibration terminée, l'écran affiche « Mesure terminée » et indique tout changement de fréquence du laser.

L'instrument peut mettre jusqu'à 12 heures pour atteindre l'équilibre thermique. Pour obtenir les meilleurs résultats, ré-exécutez l'alignement et la calibration après les 12 premières heures de fonctionnement.

- c. Exécutez le workflow Nicolet FTIR Qualification d'usine. Les workflows de vérification des performances et de qualification utilisent du polystyrène comme matériau de référence dans le spectromètre pour exécuter une série de tests standard pour vérifier le fonctionnement de votre instrument et assurer la précision de vos données.
 - i. Revenez au Tableau de bord.
 - ii. Faites défiler le panneau Workflows et sélectionnez Nicolet FTIR Qualification d'usine.
 - iii. Cliquez avec le bouton droit sur le workflow et sélectionnez **Exécuter** pour lancer le workflow.
- 6. Remettez en place tous les accessoires que vous avez retirés précédemment.

Remplacement des fenêtres du compartiment à échantillons

Le spectromètre Thermo Scientific Nicolet Summit LITE FTIR comprend un compartiment à échantillons dont les deux côtés sont dotés de fenêtres qui scellent l'instrument pour l'isoler de l'humidité et d'autres contaminants, tout en permettant au faisceau infrarouge d'entrer et de sortir du compartiment à échantillons. Les fenêtres doivent toujours être installées, même si votre système est purgé avec de l'air sec ou de l'azote.



AVERTISSEMENT





La source infrarouge située à l'intérieur de l'instrument représente une source d'inflammation. Si vous utilisez des solvants volatils, installez une hotte d'aspiration ou tout autre système de ventilation actif qui ne génère pas d'étincelles ou d'autres sources d'inflammation, et qui prévient l'accumulation de vapeurs inflammables dans l'atmosphère qui entoure l'instrument.

Nous installons des fenêtres au bromure de potassium (KBr) à l'usine, et effectuons tous nos tests de fonctionnement et de performance avec les fenêtres installées. Les fenêtres en KBr sont hygroscopiques, mais possèdent un revêtement protecteur. Sur ce spectromètre, elles fournissent une plage spectrale de travail comprise entre 8000 et 350 cm⁻¹ sans lignes d'absorption optique significatives.

Pour obtenir une performance optimale du spectromètre, les fenêtres du compartiment à échantillons doivent être transparentes (non opaques) et propres (exemptes de poussière ou d'empreintes digitales). Si vos fenêtres sont contaminées, le spectromètre peut échouer aux tests de performance et

de qualification. Vous pouvez commander de nouvelles fenêtres auprès de notre société, et les installer vous-même en suivant les instructions ci-dessous.

À NOTER

- Remplacez les fenêtres du compartiment à échantillons uniquement par des pièces de rechange que nous fournissons.
- Assurez-vous que les fenêtres n'entrent pas en contact avec des liquides.
- Les fenêtres peuvent être rayées et abîmées très facilement. Ne les touchez pas et n'essayez pas de les nettoyer. La poussière n'affectera pas le signal, tandis que les traces de doigt peuvent dégrader les performances de l'instrument et endommager de façon permanente les miroirs ou les fenêtres. Si vous souhaitez enlever la poussière à la surface d'un miroir ou d'une fenêtre, chassez-la avec un jet d'air ou d'azote propre et sec à basse pression. (N'utilisez pas d'air comprimé provenant d'un aérosol ; des contaminants risqueraient de les endommager).
- Laissez les nouvelles fenêtres dans leur emballage de protection jusqu'à ce que vous soyez prêt (e) à les insérer dans le spectromètre.
- Si vous retirez des fenêtres en KBr ou en ZnSe du spectromètre, placez-les immédiatement dans l'emballage de protection fourni (y compris le desséchant) et scellez soigneusement l'emballage.
- Les fenêtres en KBr sont transparentes ; les fenêtres en ZnSe sont jaunes.

Outils requis :

Kit de fenêtres de remplacement en ZnSe (ou en KBr) Nicolet Summit

Le kit comprend les éléments suivants :

- Fenêtres de remplacement en ZnSe (ou en KBr) (2)
- Outil de remplacement des fenêtres
- Emballage de protection (incluant le desséchant) pour le stockage des fenêtres non utilisées
- Gants en nitrile



Remplacer les fenêtres des compartiments à échantillons

1. Mise hors tension du spectromètre.

Note Si votre instrument est purgé, laissez la purge en marche pendant que vous remplacez les fenêtres pour empêcher l'air ambiant de pénétrer dans le spectromètre.

- 2. Retirez tout accessoire d'échantillonnage du spectromètre.
- 3. Retirez la première fenêtre installée.
 - a. Alignez les trois détrompeurs du bord intérieur de l'outil de remplacement des fenêtres avec les encoches du bord extérieur de la première fenêtre installée.



Note Les raccords de la fenêtre doivent être bien serrés. Stabilisez l'instrument d'une main, et retirez la fenêtre de l'autre.

b. Appuyez fermement sur l'outil et tournez-le dans le sens antihoraire pour desserrer la fenêtre.

Continuez à desserrer la fenêtre jusqu'à ce qu'elle soit libérée. Si vous inclinez l'outil vers le haut, il soutiendra la fenêtre - ce qui vous évitera de la manipuler directement.



À NOTER

Portez des gants en nitrile pour manipuler la fenêtre, et tenez-la uniquement par les bords (évitez de toucher les surfaces de la fenêtre, même si vous portez des gants).

- c. Si la fenêtre est toujours utilisable, saisissez-la par l'anneau en plastique et placez-la avec précaution dans l'emballage fourni (avec du desséchant). Rangez l'emballage dans un endroit sec et propre.
- 4. Installez la nouvelle fenêtre.
 - a. Saisissez la nouvelle fenêtre par l'anneau en plastique et placez-la avec précaution dans l'outil de remplacement de fenêtre, le pas de vis de la fenêtre tournés vers le haut.
 - b. Faites tourner la fenêtre jusqu'à ce que les trois encoches de son bord extérieur soient alignées avec les détrompeurs du bord intérieur de l'outil.
 - c. En inversant la procédure de retrait ci-dessus, inclinez l'outil et la fenêtre juste assez pour placer la fenêtre sur l'ouverture du spectromètre.
 - d. Tout en maintenant une légère pression contre la paroi du spectromètre, tournez lentement l'outil dans le sens horaire pour vous assurer que le pas de vis de la fenêtre est bien engagés.

À NOTER

La rotation de la fenêtre doit être facile au départ ; dans le cas contraire, inversez le sens de rotation et recommencez l'insertion pour éviter de fausser le pas de vis.

- e. Continuez à tourner l'outil dans le sens horaire jusqu'à ce que vous sentiez le joint torique se comprimer, puis rajouter 1/8 de tour.
- 5. Répétez les étapes 3 et 4 ci-dessus pour remplacer la deuxième fenêtre.
- 6. Mettez l'instrument sous tension, connectez-le à votre ordinateur et démarrez le logiciel OMNIC Paradigm.
- 7. Vérification des performances du spectromètre.

Avant de réaliser les étapes suivantes, assurez-vous de l'absence d'accessoires dans le compartiment échantillon.

- a. Alignez les composants optiques du spectromètre.
 - i. Dans le tableau de bord du logiciel OMNIC Paradigm, allez à Acquérir des données
 > Diagnostics > Aligner le spectromètre.
 - ii. Sélectionnez **Démarrer** pour commencer l'alignement. L'écran indique « Alignement terminé » une fois l'alignement terminé avec succès.
- b. Calibrez le laser.
 - i. Sélectionnez Acquérir des données > Diagnostic > Calibration du laser.
 - ii. Sélectionnez Démarrer pour commencer la calibration.

La calibration du laser peut prendre plusieurs minutes. Une fois la calibration terminée, l'écran affiche « Mesure terminée » et indique tout changement de fréquence du laser.

L'instrument peut mettre jusqu'à 12 heures pour atteindre l'équilibre thermique. Pour obtenir les meilleurs résultats, ré-exécutez l'alignement et la calibration après les 12 premières heures de fonctionnement.

- c. Exécutez le workflow Nicolet FTIR Qualification d'usine. Les workflows de vérification des performances et de qualification utilisent du polystyrène comme matériau de référence dans le spectromètre pour exécuter une série de tests standard pour vérifier le fonctionnement de votre instrument et assurer la précision de vos données.
 - i. Revenez au Tableau de bord.
 - ii. Faites défiler le panneau Workflows et sélectionnez Nicolet FTIR Qualification d'usine.
 - Cliquez avec le bouton droit sur le workflow et sélectionnez Exécuter pour lancer le workflow.
- 8. Si vous utilisez des tests de qualification opérationnelle (QO) pour le suivi des performances de l'instrument, nous vous recommandons de réexécuter le test de QO que vous aviez sélectionné

après avoir remplacé les fenêtres du compartiment à échantillons (en particulier, si vous avez changé le type de fenêtre). De plus amples informations sont disponibles à la rubrique « Qualification du spectromètre Summit » de l'aide en ligne.

9. Remettez en place tous les accessoires que vous avez précédemment retirés du spectromètre.