Nicolet Summit OA

Spectromètre FTIR



MANUEL DE L'UTILISATEUR

269-346500 Révision B 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 OA	
Avant la livraison du spectromètre Summit	6
Sécurité personnelle dans le cadre de l'utilisation du spectromètre Summit	
Déballage et première utilisation du spectromètre	15
Sauvegarde et récupération de vos données	
Installer un adaptateur sans fil USB	27
Comprendre la barre lumineuse du spectromètre	
Consultation de vos données via l'application OMNIC Anywhere	32

Options et accessoires	33
Installation et entretien d'un kit de purge	34

Maintenance	45
Calendrier de maintenance du spectromètre Summit	
Alignement des composants optiques du spectromètre	48
Calibration de la fréquence du laser	49
Nettoyage de l'instrument et du cristal	50
Dépose et remplacement de la plaque de cristal	
Remplacement du couvercle	54
Remplacement du desséchant	
Remplacement de la source	64

Chapitre 1

Bienvenue

Le spectromètre infrarouge à transformée de Fourier (FTIR) Thermo Scientific Nicolet Summit OA vous permet d'effectuer des analyses chimiques d'échantillons de substances en recueillant des données dans le domaine spectral de l'infrarouge moyen. 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 OA

Pour vous familiariser avec votre nouveau spectromètre, le Nicolet Summit OA, 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	
Sauvegarde et récupération de vos données	
Installer un adaptateur sans fil USB	27
Comprendre la barre lumineuse du spectromètre	
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, à l'alimentation et aux câbles de l'instrument.

- Poids de l'instrument Summit : 12,6 kg (27,8 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 OA 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.

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.

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>".



Manipulation de composants internes

purge.

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



ATTENTION



Évitez tout risque de brûlure.

provoquer un choc électrique.

Évitez tout risque de choc électrique.

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.

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

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 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.

ATTENTION



Évitez tout risque de brûlure.

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.

DANGER



Évitez tout risque d'explosion.

Avant de retirer la source du spectromètre, mettez l'instrument hors tension, débranchez le cordon d'alimentation, débranchez les lignes de purge, et retirez les accessoires et les 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.

Aperçu

Le spectromètre Nicolet Summit OA 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.

Thermo Scientific

Déballage et première utilisation du spectromètre

Le déballage et l'installation de votre spectromètre Thermo Scientific Nicolet Summit OA 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.

Le spectromètre est livré avec un couvercle anti-poussière sur le cristal ATR. Utilisez le couvercle anti-poussière lorsque vous stockez ou déplacez l'instrument, et retirez-le pour l'utilisation normale.

- Raccordez le cordon d'alimentation électrique au spectromètre et à une source de courant alternatif correctement mise à la terre.
- 3. Connectez le moniteur, le clavier et la souris.
 - a. Connectez votre moniteur au connecteur Mini DisplayPort du spectromètre.
 - b. Connectez le clavier et la souris à l'un des ports USB du spectromètre.

Note Certains dispositifs USB et Mini DisplayPort sont connus pour interférer avec le signal des dispositifs USB sans fil, tels que les claviers et les souris sans fil. Si vous remarquez des problèmes de performance au niveau de votre clavier et de votre souris sans fil (par ex., un retard ou un échec de saisie), utilisez le câble d'extension USB fourni de manière à éloigner le récepteur USB sans fil des ports USB.

4. 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.

Lorsque vous mettez l'instrument sous ou hors tension, il est possible que vous entendiez le bruit de la référence interne en polystyrène qui bouge. C'est normal.



La première fois que vous mettez le spectromètre sous tension, vous êtes invité à saisir un nouveau mot de passe. Nous vous recommandons de ne pas saisir de mot de passe à ce stade. Continuez sans saisir de mot de passe. Une fois que vous aurez terminé la configuration, vous gérerez les comptes et les mots de passe sur le dispositif en fonction des politiques de votre entreprise,

Note Si l'écran reste noir lorsque vous mettez le système sous tension, vérifiez que l'alimentation du moniteur, le connecteur DisplayPort et l'entrée du moniteur. Reportez-vous à la documentation qui accompagnait votre moniteur pour les détails de la modification des paramètres d'entrée.

5. Ouvrez le logiciel OMNIC Paradigm.

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.
 - a. Alignez les composants optiques du spectromètre.
 - i. Dans l'interface Bureau, sélectionnez 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.
 - iii. 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 Summit OA - Test PV.

Le workflow de vérification des performances utilise un matériau de référence pour exécuter une série de tests standard afin de vérifier le fonctionnement de votre instrument et d'assurer la précision de vos données.

- i. Dans le Tableau de bord de l'interface Bureau, faites défiler le panneau Workflows.
- ii. Dans le panneau Workflows, cliquez avec le bouton droit sur Nicolet Summit OA Workflow PV et sélectionnez Exécuter.

Première utilisation du spectromètre Nicolet Summit OA

~		
Name	Date Created	Last Modified
Nicolet Summit OA-PV Test	11/19/2020 1:33 PM	11/19/2020 2:14

- iii. Suivez les invites à l'écran pour terminer le workflow.
- iv. Une fois le tableau Résultats PV affiché, sélectionnez Terminer pour compléter le workflow et revenir au tableau de bord, ou sélectionnez Mesurer le spectre du squalane pour continuer avec la partie squalane en option du workflow. Mesurer un spectre de squalane fournit une référence supplémentaire pour vérifier les performances de votre système.

	thermo scientific
 Measure Squalane Spectrum Finish 	

Étapes suivantes

Félicitations ! Votre instrument est prêt à l'emploi. Cependant, avant de commencer à acquérir des données, vous devez sauvegarder votre système et créer un lecteur de récupération par mesure de précaution, afin de protéger votre système contre d'éventuels problèmes dans l'avenir. Windows 10 comprend plusieurs outils intégrés pour la sauvegarde et la récupération de vos données.

Des instructions sur la création d'un lecteur de récupération et d'une image système et sur la restauration de votre système sont disponibles à la rubrique <u>"Sauvegarde et récupération de vos</u> données".

Pour en savoir plus sur l'identification et l'analyse des échantillons avec votre spectromètre, reportezvous aux guides et tutoriels d'OMNIC Paradigm sur thermofisher.com/ftir-help.

Pour basculer sur l'interface Opérateur, allez à Visualisation/Affichage > Opérateur.

Pour les instructions détaillées de la création et de l'exécution de solutions personnalisées en utilisant l'Éditeur de Package et l'interface Opérateur, reportez-vous à « Solutions personnalisées », dans le Guide de l'utilisateur du logiciel OMNIC Paradigm.

Sauvegarde et récupération de vos données

Avant de commencer à acquérir des données, vous devez prendre quelques mesures de précaution pour protéger votre système d'éventuels problèmes futurs. Windows 10 comprend plusieurs outils intégrés pour la sauvegarde et la récupération de vos données.

IMPORTANT ! Pour vous assurer de pouvoir récupérer toutes les données en cas d'endommagement du disque dur ou du système d'exploitation, créez un support de récupération ainsi qu'une image système, et mettez en œuvre des procédures de sauvegarde de vos données.

Création d'un lecteur de récupération

Créez un lecteur de récupération de façon à pouvoir accéder aux options de démarrage avancées de Windows 10 (Récupération de l'image système, Restauration du système et Paramètres), même si votre installation de Windows est endommagée.

Matériel requis

Clé USB formatée en FAT32 avec 8 Go de stockage au minimum

À NOTER

Au cours de cette procédure, l'ensemble du contenu qui se trouve sur le disque dur sera supprimé. Avant d'exécuter cette procédure, assurez-vous d'avoir sauvegardé tous les fichiers que vous souhaitez conserver.

Créer un lecteur de réparation

- 1. Insérez la clé USB dans l'un des ports USB situés sur le côté du spectromètre.
- 2. Dans la zone de recherche de Windows 10 dans la barre des tâches, recherchez « Créer un lecteur de réparation » et sélectionnez **Créer un lecteur de réparation**.

Première utilisation du spectromètre Nicolet Summit OA



- 3. Répondez Oui à l'invite vous priant d'autoriser des modifications sur votre équipement.
- 4. Désélectionnez « Sauvegarder les fichiers système sur le lecteur de récupération » et cliquez sur Suivant.
- 5. Sélectionnez votre lecteur dans la liste des lecteurs disponibles et cliquez sur Suivant.



- 6. Cliquez sur Créer. Cliquez sur Terminer quand la tâche est terminée afin de fermer la fenêtre.
- 7. Retirez la clé USB et stockez-la en lieu sûr. Nous vous conseillons d'étiqueter la clé « Lecteur de récupération Summit ».

Pour en savoir plus sur la création de supports de récupération, consultez l'aide en ligne de Microsoft.

Création d'une image système

Une image système est une capture instantanée de l'ensemble de votre système - y compris vos données et fichiers, applications installées et autres paramètres. En cas de défaillance complète du disque dur ou du système d'exploitation, vous pouvez utiliser l'image système pour restaurer votre dispositif à l'état dans lequel il se trouvait lorsque l'image a été créée. C'est à vous de décider de la

fréquence de mise à jour de l'image système, mais il est conseillé de conserver une copie propre de l'image à l'installation.

En fonction de la quantité de données que vous copierez dans l'image système, le processus peut prendre un certain temps. La copie d'une large quantité de données peut prendre plus d'une heure.

Matériel requis

• SSD au format NTFS

Créer une image système

- 1. Raccordez votre SSD à l'un des ports USB du spectromètre.
- Dans Windows 10, ouvrez le Panneau de configuration et sélectionnez Sauvegarder et restaurer (Windows 7).
- 3. Dans le panneau de gauche, sélectionnez Créer une image système.

🛞 Backup and Restore (Windows 7)				
← → ✓ ↑ 🐌 > Control Panel > System and Security > Backup and Restore (Windows 7)				
Control Panel Home	Back up or restore your files			
🗣 Create a system image	Backup			
💡 Create a system repair	disc Windows Backup has not been set up.	Set up backup		
	Restore Windows could not find a backup for this computer.			

4. Sélectionnez Sur un disque dur et choisissez votre SSD dans la liste. Cliquez sur Suivant.

Première utilisation du spectromètre Nicolet Summit OA

👘 Create a system image	
Where do you want to save the backup?	
A system image is a copy of the drives required for Windows to run. It can also include additional drives. A system image can be used to restore your computer if your hard drive or computer ever stops working: however, you can't choose individual items to restore.	
On a hard disk	
🔜 Backup (E:) 465.60 GB free 🗸 🗸	
O On one or more DVDs	
O on one or more DVDs	
O On one or more DVDs	
On one or more DVDs	
On one or more DVDs DVD Drive (D:) On a network location	
On one or more DVDs DVD Drive (D:) On a network location Select	
On one or more DVDs DVD Drive (D:) On a network location Select	
On one or more DVDs DVD Drive (0:) On a network location Select	
On one or more DVDs DVD Drive (D:) On a network location Select	
On one or more DVDs DVD Drive (D:) On a network location Select	

- 5. Confirmez les paramètres de sauvegarde et cliquez sur Démarrer la sauvegarde.
- 6. À l'invite « Voulez-vous créer un disque de réparation système ? », cliquez sur Non.
- 7. Cliquez sur **Fermer** et retirez le SSD du Summit en toute sécurité. Étiquetez le lecteur « Sauvegarde du Summit », et stockez-le en lieu sûr.

Restaurer votre système

En cas de défaillance du système d'exploitation, restaurez votre système à l'aide de votre lecteur de récupération et de l'image système.

Matériel requis

- Lecteur de récupération USB
- SSD avec votre image système
- Restaurer votre système à l'aide du lecteur de récupération et de l'image système
- 1. Éteignez le spectromètre Summit.
- 2. Raccordez un clavier externe à l'un des ports USB du spectromètre Summit.
- 3. Raccordez le lecteur de récupération USB à l'un des ports USB.
- 4. Rallumez le système.
- 5. Lorsque l'écran de démarrage Thermo Scientific s'affiche, appuyez plusieurs fois sur la touche

Echap du clavier pour ouvrir le BIOS du système.

Si une boîte de dialogue affiche le message « Quit Without Saving ? », sélectionnez No à l'aide des flèches du clavier.

- 6. Accédez à l'onglet Save & Exit à l'aide des flèches du clavier.
- 7. Naviguez à travers la section Boot Override et sélectionnez votre clé USB de récupération. Appuyez sur la touche **Enter** pour démarrer à partir de l'image de récupération.



- 8. Lorsque l'écran « Choisir la disposition du clavier » apparaît, retirez la clé USB de récupération du port USB, et insérez le SSD avec l'image système.
- 9. Sélectionnez la disposition du clavier en fonction de votre préférence.
- 10. Dans l'écran Choisir une option, sélectionnez Dépannage.



- 11. Sélectionnez Options avancées.
- 12. Dans Options avancées, sélectionnez Récupération de l'image système.

Première utilisation du spectromètre Nicolet Summit OA



- 13. Sélectionnez Windows 10 comme système d'exploitation cible.
- 14. Dans la fenêtre « Sélectionner une sauvegarde d'image système », utilisez la dernière image système disponible ou sélectionnez une autre image système. Cliquez sur **Suivant**.



15. Cliquez sur **Suivant** pour conserver l'état par défaut des options de restauration supplémentaires, puis sur **Terminer** pour restaurer l'image système. Cliquez sur **Oui** pour confirmer.

La restauration prend environ 15 minutes et le système redémarrera automatiquement une fois l'opération terminée.

Une fois le système restauré, retirez le SSD du spectromètre en toute sécurité, et stockez le lecteur de récupération et le SSD de l'image système dans un endroit sûr.

De plus amples informations sur la restauration du système à partir de l'image système sont disponibles dans l'aide en ligne de Microsoft.

Sauvegarde des données de votre système

Outre la création de supports de récupération et d'une image système, vous devez régulièrement effectuer une sauvegarde de vos données. Windows 10 propose plusieurs options de sauvegarde des

Première utilisation du spectromètre Nicolet Summit OA

fichiers, dont l'outil Historique des fichiers. Les informaticiens de votre société doivent déterminer les méthodes de sauvegarde les mieux adaptées à vos besoins.

Installer un adaptateur sans fil USB

Vous pouvez connecter votre spectromètre FTIR Nicolet Summit OA à un réseau sans fil en utilisant l'adaptateur sans fil USB en option. Si vous l'avez acheté avec votre spectromètre, l'adaptateur est livré préinstallé.

Pour installer l'adaptateur sans fil USB, insérez l'adaptateur dans un port USB. Pour une installation plus permanente, installez l'adaptateur dans un port sous le couvercle de l'instrument.

Installer un adaptateur sans fil USB

1. Insérez l'adaptateur dans un port USB disponible.



2. Le système reconnaîtra automatiquement l'adaptateur. Pour les informations pour la connexion à un réseau en utilisant votre adaptateur, consultez l'assistance de Microsoft.

Installer le système en utilisant un port USB sous le couvercle de l'instrument

Pour ce type d'installation, vous devez retirer la plaque du cristal et le couvercle de l'instrument.

- 1. Pour retirer la plaque du cristal et le couvercle de l'instrument, suivez les instructions de la rubrique "Remplacement du couvercle"
- 2. Le couvercle déposé, insérez l'adaptateur dans l'un ou l'autre des ports USB disponibles.



3. Remettez en place le couvercle de l'instrument et la plaque du cristal.

Votre instrument est prêt à être connecté à un réseau sans fil.

Comprendre la barre lumineuse du spectromètre

Le spectromètre Thermo Scientific Nicolet Summit OA 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 OA



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 2-1 : OMNIC Anywhere dans le navigateur web



Un spectromètre Nicolet Summit OA 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> ».

Options et accessoires

Le Nicolet Summit OA peut être utilisé avec le kit de purge en option.

stallation et entretien d'un kit de purge

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

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

- Une clé anglaise de 3/4"
- Une clé anglaise de 11/16"
- Un tournevis cruciforme n° 2
- Clé hexagonale de 3/32"

Options et accessoires

- Clé hexagonale de 7/64"
- 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", 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" 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.
- 2. Mettez hors tension le spectromètre.
 - a. Pour éteindre le spectromètre, appuyez sur le bouton d'alimentation.
 - b. Débranchez le cordon d'alimentation.
- 3. "Retrait de la plaque de cristal"
- 4. Déposez le couvercle de l'instrument.
 - a. Inclinez délicatement l'instrument sur son dos pour qu'il repose sur les ailettes verticales du dissipateur thermique.
 - b. En utilisant une clé de 7/64", desserrez les quatre vis qui fixent le couvercle à la base de l'instrument.



- c. Inclinez délicatement l'instrument pour le remettre sur ses pieds. Lorsque vous le remettez à la verticale, les quatre vis retombent. Faites attention à ne pas les perdre.
- d. Soulevez droit le couvercle de la base.



5. Retirez le support de desséchant.

Options et accessoires

a. 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.



b. Retirez et jetez le grand joint torique.



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

- 6. Ouvrez le kit de purge et le sachet de desséchant scellé.
- 7. 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.

8. 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.



- 9. Faites passer la ligne de purge à travers l'ouverture d'entrée de purge à l'arrière de l'instrument.
 - a. Abaissez délicatement le couvercle sur la base tout en faisant passer le tube de purge dans l'encoche du couvercle.



- b. Abaissez complètement le couvercle. Lorsque vous abaissez le couvercle, le tube de purge est bloqué dans l'encoche entre le couvercle et le dissipateur thermique.
- c. Le tube de purge étant repoussé sur le côté, inclinez l'instrument sur le dos pour qu'il repose sur les ailette du dissipateur thermique.
- d. Insérez les quatre vis retirées au préalable. Serrez les quatre vis.
- e. Inclinez l'instrument pour le remettre sur ses pieds.
- f. Faites entrer la tubulure en excès à l'intérieur l'instrument, ou sortez-la en si nécessaire.



10. "Remplacement de la plaque en cristal".

- 11. Raccordez le coupleur de débit de l'ensemble de tuyauterie murale (étiqueté « Vers le spectromètre ») à la tubulure de purge.
- 12. 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).

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

- 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.



- 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).
- 3. Réglez le débitmètre du spectromètre sur 1 scfh (0,47 l/min).

À NOTER

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

4. 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.

[Cette page est intentionnellement laissée vide]

Chapitre 4

Maintenance

Votre spectromètre Nicolet Summit OA 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	
Alignement des composants optiques du spectromètre	
Calibration de la fréquence du laser	
Nettoyage de l'instrument et du cristal	
Dépose et remplacement de la plaque de cristal	
Remplacement du couvercle	54
Remplacement du desséchant	58
Remplacement de la source	64

Calendrier de maintenance du spectromètre Summit

Votre spectromètre Thermo Scientific Nicolet Summit OA 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 Paradigm comprend un workflow de vérification des performances conçu pour votre spectromètre. 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.

Nous vous recommandons d'exécuter le workflow de vérification des performances (PV) 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

Maintenance

desséchant qui absorbent l'humidité. Surveillez l'humidité de l'instrument dans le logiciel OMNIC Paradigm.

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 4-1 : 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 du spectromètre si le workflow Nicolet Summit OA - Test PV indique que le test a échoué. L'alignement du spectromètre optimise l'énergie qui atteint le détecteur et maximise le signal du détecteur. Si un test de performance 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 [100] pour ouvrir la vue 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

Calibrez la fréquence de votre laser si le workflow Nicolet Summit OA - Test PV 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 de l'instrument et du cristal

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

Nettoyage du spectromètre

Nettoyez délicatement l'extérieur de l'instrument avec un chiffon propre et doux légèrement humidifié avec un solvant doux, par exemple du Windex. Vaporisez le produit sur le chiffon et ne nettoyez que la surface extérieure peinte. Veillez à ne pas répandre de solvant sur les composants électroniques à l'arrière de l'instrument.

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.

Nettoyage de la plaque en cristal ATR

Utilisez un chiffon ou un essuie-tout propre et sec non-abrasif pour nettoyer le cristal ATR.

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.

Enfilez des gants de protection pour nettoyer ou changer la plaque en cristal.

- 1. En utilisant un chiffon ou un essuie-tout en papier propre, absorbez le liquide s'il y en a sur le cristal. Jetez l'essuie-tout en papier usagé.
- 2. Avec une pipette, appliquez quelques gouttes d'heptane sur le cristal et tout autour de celui-ci.
- 3. Avec un nouvel essuie-tout en papier propre, essuyez le cristal sur toute sa longueur pour éliminer tout liquide éventuel. Essuyez bien la zone autour du cristal.



4. Répétez si nécessaire et finissez toujours avec un chiffon ou un essuie-tout en papier propre et sec.

Dépose et remplacement de la plaque de cristal

La plaque de cristal doit être changée tous les 12 à 18 mois ou selon les besoins.

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.

Enfilez des gants de protection pour nettoyer ou changer la plaque en cristal.

À NOTER

Veillez à ce que rien ne tombe de l'instrument quand le couvercle est déposé.

MATÉRIEL REQUIS

Clé hexagonale de 3/32"

Retrait de la plaque de cristal

- 1. Retirez le couvercle anti-poussière du cristal.
- 2. En utilisant la clé hexagonale de 3/32", desserrez complètement les trois vis qui fixent la plaque en cristal à l'instrument.

Maintenance



- 3. Soulevez droit la plaque en cristal de l'instrument.
- 4. Utilisez un chiffon propre et sec pour nettoyer le pourtour de la bride, sous la zone qui accueille la plaque.

Remplacement de la plaque en cristal

- 1. Mettez la nouvelle plaque en place sur la bride.
- 2. Placez les rondelles sur les vis et insérez les vis dans les trois trous. Les vis doivent descendre droit dans les trous.
- 3. Tout en appuyant sur la plaque en cristal, serrez les deux vis sur les deux côtés de la plaque, puis serrez la troisième vis (la plus proche du dos de l'instrument). Serrez les vis manuellement, mais à fond.

Si une vis tourne librement, faites-la remonter puis réinsérez-la pour corriger l'alignement.

Remplacement du couvercle

OUTILS REQUIS

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

- Une clé hexagonale de 7/64"
- Une clé hexagonale de 3/32"

Remplacer le couvercle du spectromètre

- 1. Mise hors tension du spectromètre.
 - a. Pour éteindre le spectromètre, appuyez sur le bouton d'alimentation.
 - b. Débranchez le cordon d'alimentation.
 - c. Arrêtez la purge, si applicable, et débranchez la ligne de purge de l'instrument.
- 2. Retirez la plaque en cristal.
 - a. Retirez le couvercle anti-poussière du cristal.
 - b. Desserrez complètement les trois vis qui fixent la plaque en cristal à l'instrument.



- c. Soulevez droit la plaque en cristal de l'instrument.
- d. Utilisez un chiffon propre et sec pour nettoyer le pourtour de la bride, sous la zone qui accueille la plaque.
- 3. Déposez le couvercle de l'instrument.
 - a. Inclinez délicatement l'instrument sur son dos pour qu'il repose sur les ailettes verticales du dissipateur thermique.
 - b. En utilisant une clé de 7/64", desserrez les quatre vis qui fixent le couvercle à la base de l'instrument.



c. Inclinez délicatement l'instrument pour le remettre sur ses pieds. Lorsque vous le remettez à la verticale, les quatre vis retombent. Faites attention à ne pas les perdre.

d. Soulevez droit le couvercle de la base.



- 4. Remettez en place le couvercle de l'instrument.
 - a. Posez délicatement le couvercle sur la base de l'instrument.
 - b. Inclinez l'instrument sur le dos pour qu'il repose sur les ailette du dissipateur thermique.
 - c. Insérez les quatre vis retirées au préalable. Serrez toutes les vis.
 - d. Inclinez l'instrument pour le remettre sur ses pieds.
- 5. Remplacez la plaque en cristal.
 - a. Remettez la plaque en place sur la bride.
 - b. Placez les rondelles sur les vis et insérez les vis dans les trois trous. Les vis doivent descendre droit dans les trous.
 - c. Tout en appuyant sur la plaque en cristal, serrez les deux vis sur les deux côtés de la plaque, puis serrez la troisième vis (la plus proche du dos de l'instrument). Serrez les vis manuellement, mais à fond.

Si une vis tourne librement, faites-la remonter puis réinsérez-la pour corriger l'alignement.

6. Mise sous tension du spectromètre.

Maintenance

- 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.
- 7. Vérification des performances du spectromètre.
 - 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. Calibrer 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 Summit OA Test PV**. Le workflow de vérification des performances utilise un matériau de référence 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 Summit OA Test PV.
 - iii. Cliquez avec le bouton droit sur le workflow et sélectionnez **Exécuter** pour lancer le workflow.

Remplacement du desséchant

Surveillez l'humidité de l'instrument en utilisant le logiciel OMNIC Paradigm. Changez le desséchant comme requis ou en fonction des politiques de votre entreprise.

Pour remplacer le desséchant, vous devez retirer le couvercle de l'instrument, retirer le support de desséchant et remplacer les cartouches ainsi que le joint torique.

OUTILS REQUIS

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

- Une clé hexagonale de 7/64"
- Un kit de remplacement du desséchant
- Une clé hexagonale de 3/32"
- 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 de déposer le couvercle de l'instrument et d'ouvrir le compartiment à desséchant, mettez l'instrument hors tension, débranchez le cordon d'alimentation, et enlevez tous 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. Arrêtez la purge, si applicable, et débranchez la ligne de purge de l'instrument.

Note Vous pouvez utiliser un support de desséchant avec ou sans connecteur de purge pour maintenir l'humidité requise. Les illustrations ci-dessous présentent une version de support de desséchant qui n'a pas été purgée. Les instructions pour le remplacement des cartouches de desséchant sont identiques pour les deux types de support.

- 2. <u>"Retrait de la plaque de cristal"</u>
- 3. Déposez le couvercle de l'instrument.
 - a. Inclinez délicatement l'instrument sur son dos pour qu'il repose sur les ailettes verticales du dissipateur thermique.
 - b. En utilisant une clé de 7/64", desserrez les quatre vis qui fixent le couvercle à la base de l'instrument.



c. Inclinez délicatement l'instrument pour le remettre sur ses pieds. Lorsque vous le remettez à la verticale, les quatre vis retombent. Faites attention à ne pas les perdre.

d. Soulevez droit le couvercle de la base.



- 4. Retirez le support de desséchant.
 - a. 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.



b. Retirez et jetez le grand joint torique.

Maintenance



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

5. Retirez les cartouches de desséchant et remplacez l'indicateur d'humidité.



Le contenu des cartouches de desséchant (tamis moléculaire d'aluminosilicate métallique hydraté) peut être nocif s'il est ingéré. Si vous jetez les réservoirs de desséchant saturés, veillez à les éliminer de manière appropriée.

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.



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.

- 6. 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.
- 7. 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.

- 8. Remettez en place le couvercle de l'instrument.
 - a. Posez délicatement le couvercle sur la base de l'instrument.
 - b. Inclinez l'instrument sur le dos pour qu'il repose sur les ailette du dissipateur thermique.
 - c. Insérez les quatre vis retirées au préalable. Serrez toutes les vis.

Maintenance

- d. Inclinez l'instrument pour le remettre sur ses pieds.
- e. "Remplacement de la plaque en cristal".
- 9. 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.
- 10. Vérification des performances du spectromètre.
 - 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. Calibrer 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 Summit OA Test PV**. Le workflow de vérification des performances utilise un matériau de référence 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 Summit OA Test PV.
 - 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 a été purgé, débranchez les lignes de purge du spectromètre. De plus amples informations sont disponibles à la rubrique <u>"Installation et entretien d'un kit de purge"</u>.

ATTENTION



d.

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



b. 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.
- 5. Vérification des performances du spectromètre.
 - 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. Calibrer 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 Summit OA - Test PV**. Le workflow de vérification des performances utilise un matériau de référence pour exécuter une série de tests standard pour vérifier le fonctionnement de votre instrument et assurer la précision de vos données.

Maintenance

- i. Revenez au Tableau de bord.
- ii. Faites défiler le panneau Workflows et sélectionnez Nicolet Summit OA Test PV.
- iii. Cliquez avec le bouton droit sur le workflow et sélectionnez **Exécuter** pour lancer le workflow.

[Cette page est intentionnellement laissée vide]