Nicolet Summit OA

Espectrómetro FTIR



GUÍA DEL USUARIO

269-346500 Revisión B Abril de 2021

© 2020 Thermo Fisher Scientific Inc. Reservados todos los derechos.

Para obtener soporte técnico, póngase en contacto con: www.thermofisher.com

Thermo Fisher Scientific Inc. suministra este documento a sus clientes con la compra de un producto para que lo apliquen a su funcionamiento. Este documento está protegido por derechos de autor y su reproducción total o parcial está estrictamente prohibida, salvo con la autorización por escrito de Thermo Fisher Scientific Inc.

El contenido de este documento está sujeto a cambios sin previo aviso. Toda la información técnica de este documento se incluye solo para consulta. Las especificaciones y configuraciones del sistema descritas en este documento sustituyen a toda la información anterior que haya recibido el comprador.

Thermo Fisher Scientific Inc. no se manifiesta sobre la exactitud, precisión o ausencia de errores del presente documento ni será responsable de posibles errores, omisiones, daños o pérdidas que puedan derivarse de su empleo, aun cuando la información contenida en él se siga de forma correcta.

Este documento no forma parte de ningún contrato de venta entre Thermo Fisher Scientific Inc. y el comprador. Este documento no regirá ni modificará en ningún caso los Términos y condiciones de venta, que prevalecerán en caso de conflicto entre la información de ambos documentos.

Solo para uso en investigación. Este instrumento o accesorio no es un producto sanitario y no está diseñado para utilizarse en la prevención, el diagnóstico, el tratamiento o la cura de enfermedades.

ADVERTENCIA



Evite el peligro de fuego o explosiones.

Este instrumento o accesorio no está diseñado para utilizarse en ambientes explosivos.

Índice

| Bienvenida | 1 |
|----------------------------------|---|
| Símbolos informativos utilizados | 2 |
| Pedir componentes | 3 |
| Contactar con nosotros | 3 |

| Primeros pasos con el espectrómetro Nicolet Summit OA | 5 |
|---|----|
| Antes de recibir el espectrómetro Summit | 6 |
| Seguridad personal con el espectrómetro Summit | 9 |
| Desembalar y preparar el espectrómetro | 15 |
| Copia de seguridad y recuperación de datos | 20 |
| Instalar un adaptador inalámbrico USB | |
| Funcionamiento de la barra luminosa del espectrómetro | |
| Ver los datos con la aplicación OMNIC Anywhere | |
| | |

| Opciones y accesorios | |
|-------------------------------------|----|
| Instalar y mantener un kit de purga | 34 |

| Mantenimiento | 45 |
|--|----|
| Programa de mantenimiento del espectrómetro Summit | |
| Alinear la óptica del espectrómetro | 48 |
| Calibrar la frecuencia del láser | 49 |
| Limpiar el instrumento y el cristal | 50 |
| Quitar y sustituir la placa de cristal | 52 |
| Sustituir la cubierta | 54 |
| Sustituir el desecante | 58 |
| Sustituir la fuente | 64 |

Capítulo 1

Bienvenida

El espectrómetro de infrarrojo por transformada de OA Fourier (FTIR) Nicolet Summit OA de Thermo Scientific permite realizar análisis químicos de muestras de materiales mediante la adquisición de datos en el rango espectral del IR medio. El sistema integra funciones de verificación, un potente paquete de software y muchas otras características que facilitan la adquisición de datos. La instalación de hardware opcional y los procedimientos de mantenimiento y reparación puede realizarlos el usuario. En este documento, u otra documentación provista, encontrará toda la información necesaria.

Asegúrese de leer la guía de seguridad que acompaña al sistema antes de utilizar el instrumento.

Nota Se recomienda mantener el instrumento sellado y desecado o purgado en todo momento. Los daños causados al equipo por no mantener sellado y desecado o purgado el instrumento no están cubiertos por la garantía. Si tiene alguna duda a este respecto, póngase en contacto con nosotros.

Símbolos informativos utilizados

Las precauciones de seguridad y otra información importante se presentan en el siguiente formato:

PELIGRO



Peligro.Indica una situación peligrosa que, de no evitarse, ocasionará lesiones graves o mortales.

ADVERTENCIA



Peligro. Indica una situación peligrosa que, de no evitarse, puede ocasionar lesiones graves o mortales.

PRECAUCIÓN



Peligro.Indica una situación peligrosa que, de no evitarse, puede ocasionar lesiones leves o moderadas.

AVISO

Siga las instrucciones marcadas con esta etiqueta para evitar que se produzcan daños en el hardware del sistema o pérdida de datos.

Nota Contiene información complementaria muy útil.

Bienvenida

Pedir componentes

Para pedir componentes, póngase en contacto con nosotros.

Si necesita enviarnos el instrumento o un accesorio para su reparación, llámenos o envíenos un correo electrónico primero para conocer los requisitos del envío y otras instrucciones.

Contactar con nosotros

Para obtener Soporte Técnico, póngase en contacto con: www.thermofisher.com

[Página en blanco]

Primeros pasos con el espectrómetro Nicolet Summit OA

Para poder trabajar con el nuevo espectrómetro Nicolet Summit OA, es preciso revisar los requisitos de seguridad y emplazamiento antes de que llegue el instrumento, de desembalarlo y prepararlo, y de conectarle los accesorios adicionales.

| Antes de recibir el espectrómetro Summit | 6 |
|---|----|
| Seguridad personal con el espectrómetro Summit | 9 |
| Desembalar y preparar el espectrómetro | 15 |
| Copia de seguridad y recuperación de datos | |
| Instalar un adaptador inalámbrico USB | |
| Funcionamiento de la barra luminosa del espectrómetro | |
| Ver los datos con la aplicación OMNIC Anywhere | 31 |

Antes de recibir el espectrómetro Summit

Antes de recibir su espectrómetro, prepare un área de trabajo idónea para garantizar la mayor exactitud de los datos y el mejor rendimiento a largo plazo.

En esta sección se explica cómo proceder a la llegada del espectrómetro y cómo los factores ambientales y eléctricos afectan al rendimiento del espectrómetro. Para obtener una explicación más exhaustiva sobre los requisitos, consulte la Información sobre seguridad y emplazamiento del espectrómetro.

Desembalar el espectrómetro

Antes de abrir la caja, debe seguir dos pasos importantes al recibir el instrumento:

• Revisar el exterior de la caja de transporte por si está dañada

Si observa daños externos, póngase en contacto con nosotros o con el distribuidor local para solicitar instrucciones.

• Dejar que el espectrómetro alcance la temperatura ambiente

Dentro de la caja de transporte, el espectrómetro se encuentra en una bolsa de plástico cerrada herméticamente que lo mantiene seco. Deje pasar 24 horas antes de abrir la bolsa, para que el espectrómetro alcance la temperatura ambiente de la sala. Si abre la bolsa antes de que el espectrómetro se haya calentado, podría condensarse humedad en los componentes ópticos y dañarlos de forma irreversible.

La garantía no cubre daños causados por métodos de desplazamiento inadecuados o por retirar la bolsa de plástico hermética antes de que el instrumento haya alcanzado la temperatura ambiente.

Preparar el área de trabajo

Antes de que llegue el instrumento, asegúrese de que el área de trabajo soporte con seguridad el espectrómetro. Además del espacio necesario para colocar el espectrómetro, tenga en cuenta los distintos requisitos ambientales y eléctricos.

Dimensiones del espectrómetro

El espectrómetro Summit necesita muy poco espacio, pero se debe dejar sitio a su alrededor para que el calor de los orificios de ventilación se pueda disipar y sea fácil acceder al interruptor de encendido/apagado del instrumento y a sus puertos y cables.

- Peso del instrumento Summit: 12,6 kg (27,8 lb)
- Dimensiones (An x Al x Pr): 34 x 24 x 32 cm (13,3 x 9,6 x 12,7 pulg.)

Factores ambientales

El espectrómetro Summit es un instrumento resistente diseñado para utilizarse en muchos entornos. No obstante, para optimizar su rendimiento, conviene mantenerlo en ambientes con poca humedad y escaso polvo. El espectrómetro funciona con fiabilidad a temperaturas de entre 15 y 35 °C, pero su rendimiento óptimo se asegura a temperaturas de entre 20 y 22 °C.

La humedad puede condensarse en el interior del instrumento y deteriorar sus componentes internos. Algunas medidas de precaución ayudan a proteger el instrumento de la humedad:

- Conservar el desecante del instrumento, incluso cuando esté almacenado.
- Evitar cambios bruscos de temperatura.
 - Mantener el instrumento alejado de fuentes de aire frío o caliente, como aparatos de calefacción o aire acondicionado o grandes ventanales.

Si el espectrómetro va a estar ubicado en un ambiente especialmente húmedo, considere instalar un kit de gas de purga.

Requisitos eléctricos

La alimentación eléctrica del espectrómetro debe proceder de fuentes de alimentación dedicadas, ininterrumpidas y libres de lo siguiente:

- Caídas de tensión
- Picos de corriente
- Cambios de frecuencia
- Otras perturbaciones eléctricas

Si sospecha que hay problemas eléctricos, recomendamos realizar una inspección de la calidad de la alimentación eléctrica. Póngase en contacto con nosotros o con la compañía eléctrica local para obtener más información.

Especificaciones del suministro eléctrico

En la siguiente tabla se muestran las especificaciones del servicio eléctrico. Póngase en contacto con nuestro representante de servicio de su zona si tiene alguna pregunta sobre los requisitos.

| Requisitos | Especificación |
|----------------------------|---|
| Corriente de entrada | 1,5 A máx. |
| Tensión de entrada | De 100 a 240 V CA |
| Frecuencia de línea | De 50 a 60 |
| Perturbaciones de línea | Las flechas de tensión, sobretensiones y otras perturbaciones eléctricas no deben superar el 10 % del valor de la tensión de entrada. |
| Ruido | Menos de 2 V (modo común) Menos de 20 V (modo normal) |

Seguridad personal con el espectrómetro Summit

Aunque el diseño del espectrómetro Nicolet Summit OA lo hace un instrumento seguro, debe tomar ciertas medidas de precaución para protegerse de peligros potenciales que pueden sufrirse durante su uso y mantenimiento habituales.

PRECAUCIÓN



En esta guía se describen algunos peligros potenciales que deben tenerse en cuenta, pero no es una guía exhaustiva. Antes de utilizar el instrumento, consulte la Información sobre seguridad y emplazamiento para obtener una descripción más completa de los peligros potenciales.

Peligros potenciales durante el uso normal

Durante el uso normal, los peligros proceden en su mayoría de:

- Muestras y disolventes potencialmente peligrosos
- Exposición a seleniuro de zinc (ZnSe)
- Piezas del espectrómetro calientes o emisoras de calor
- Láser del instrumento

Para intentar evitar lesiones personales o daños al instrumento, debe conocer los posibles peligros y tomar algunas precauciones.

Muestras y disolventes peligrosos

Tome medidas de precaución especiales si utiliza o planea medir muestras o disolventes potencialmente peligrosos, como gases presurizados o disolventes corrosivos o inflamables.

USAR VENTILACIÓN SUFICIENTE

El espectrómetro no tiene requisitos especiales de ventilación, pero con ciertos tipos de análisis puede ser necesario aumentar la ventilación. Asegúrese de contar con la ventilación necesaria cuando vaya a analizar muestras altamente tóxicas, disolver muestras en disolventes que pueden interaccionar con la fuente de infrarrojos, o medir gases inflamables. La pirólisis de los disolventes que contienen hidrocarburos halogenados puede producir ácido clorhídrico (HCI), ácido fluorhídrico (HF) o fosgeno (COCI₂).

ADVERTENCIA



Evite la inhalación de sustancias tóxicas. El ácido clorhídrico, el ácido fluorhídrico y el fosgeno son altamente tóxicos. Si utiliza disolventes que contienen hidrocarburos halogenados, asegúrese de que el área de trabajo esté debidamente ventilada.

DISOLVENTES VOLÁTILES E INFLAMABLES

La fuente de infrarrojos del interior del espectrómetro puede prender las muestras y disolventes volátiles e inflamables. Cuando trabaje con disolventes y muestras inflamables, siga estas indicaciones:

- Trabaje con las ventanas del compartimento de muestras instaladas.
- Asegúrese de que el área de trabajo esté debidamente ventilada y cuente con un sistema de ventilación activo que no permita chispas ni otras fuentes de ignición y que impida la acumulación de vapores inflamables en los alrededores del instrumento.
- No deje muestras ni disolventes inflamables cerca del instrumento.
- No deje muestras ni disolventes inflamables en el compartimento de muestras más tiempo del necesario.
- Purgue el espectrómetro con nitrógeno o aire seco limpio.

DISOLVENTES CORROSIVOS

Los disolventes que producen vapores de HCI o HF en el compartimento de muestras pueden dañar seriamente el sistema. Si utiliza disolventes halogenados, purgue el instrumento con nitrógeno o aire seco limpio.

AVISO

La garantía no cubre daños del equipo causados por falta de purga del instrumento.

PELIGRO BIOLÓGICO O MATERIALES RADIACTIVOS Y AGENTES INFECCIOSOS

Las muestras biológicas, tales como tejidos, líquidos corporales, agentes infecciosos y sangre, pueden transmitir enfermedades infecciosas. Siga los protocolos del Programa de bioseguridad de su organización para el trabajo con materiales potencialmente infecciosos.

Materiales tóxicos

El espectrómetro Nicolet Summit OA incluye un cristal ATR de seleniuro de zinc (ZnSe); otros instrumentos Summit podrían utilizar ventanas del compartimento de muestras opcional de ZnSe.

ADVERTENCIA



Evite la inhalación e ingestión de sustancias tóxicas.

El seleniuro de zinc (ZnSe) es tóxico. Consulte las medidas de control de manipulación y exposición en la Ficha de datos de seguridad de ZnSe del fabricante, en www.specac.com.

Fuentes de calor

Los componentes del espectrómetro pueden alcanzar temperaturas muy altas durante el funcionamiento normal. Tenga especial cuidado en torno a la fuente de infrarrojos y los orificios de ventilación del espectrómetro.

La superficie de la fuente de infrarrojos que da al exterior se encuentra en la base del instrumento y puede alcanzar temperaturas elevadas. No toque la parte inferior del instrumento durante el funcionamiento ni durante un rato después de utilizarlo.

La salida de ventilación del espectrómetro se encuentra en el panel izquierdo del instrumento. Durante el uso normal, el espectrómetro disipa el aire caliente de sus componentes internos hacia el exterior del espectrómetro a través de estos orificios de ventilación. Deje espacio suficiente alrededor de los orificios de ventilación del instrumento para que se pueda disipar el aire caliente.

Seguridad del láser y óptica

Durante el uso normal del espectrómetro, el usuario no se encuentra nunca expuesto a niveles de radiación láser inseguros. Si se retira la cubierta durante un procedimiento de servicio, podría necesitar tomar precauciones especiales, como utilizar gafas de protección. Si son necesarias, el personal técnico le informará.

ADVERTENCIA



Evite lesiones personales.

Jamás mire directamente al rayo láser ni a su reflejo. No modifique nunca el láser, incluso si debe sustituir un láser defectuoso; podría quedar expuesto a la luz láser o a altos voltajes.

Peligros potenciales durante el mantenimiento

Los peligros a los que puede estar expuesto durante las tareas de mantenimiento del instrumento son distintos a los del uso normal del mismo. Durante el mantenimiento, los principales peligros se encuentran en las tareas de purga del instrumento y de manipulación de sus componentes internos.

Purga del instrumento

En ambientes especialmente húmedos, recomendamos instalar una fuente de nitrógeno o aire seco limpio para purgar el espectrómetro. Purgar el instrumento ayuda a proteger los componentes ópticos internos contra daños derivados por el exceso de humedad y de disolventes corrosivos, además de asegurar resultados más exactos.

Consulte <u>"Instalar y mantener un kit de purga"</u> para conocer más detalles sobre la adquisición e instalación de un kit de purga para el espectrómetro.

PELIGRO

Evite el peligro de fuego y explosiones.



- Utilice solo nitrógeno o aire seco para purgar el espectrómetro.
- No utilice nunca un gas tóxico, inflamable o combustible para purgar el instrumento. El gas de purga debe estar libre de aceite y otros materiales reactivos. El calor producido por la fuente o por la absorción de láser puede encender los gases inflamables o los materiales reactivos que se encuentran en el gas de purga.

Trabajo con componentes internos

Por lo general, no debería necesitar retirar la cubierta del instrumento ni intentar sustituir componentes internos. No obstante, en caso de que necesite retirar la cubierta para alguna tarea de mantenimiento, recuerde que corre el riesgo de descargas, quemaduras y exposición a la luz láser.

PRECAUCIÓN



Evite el peligro de descargas eléctricas.

Incluso después de haberlo desconectado de todas las fuentes de tensión, los condensadores del instrumento pueden permanecer cargados durante un máximo de 30 segundos y pueden provocar una descarga eléctrica.

PRECAUCIÓN



Evite el peligro de quemaduras.

Los componentes internos, en particular la fuente de infrarrojos, pueden alcanzar temperaturas extremadamente altas durante el funcionamiento normal. Apague el instrumento y espere al menos 10 minutos antes de sustituir componentes.

ADVERTENCIA



Evite lesiones personales.

- Jamás mire directamente al rayo láser ni a su reflejo. No modifique nunca el láser. Podría quedar expuesto a luz láser o a altos voltajes.
- Si ajusta el láser o realiza procedimientos no descritos en las guías o manuales del usuario, podría quedar expuesto a radiaciones peligrosas.

SUSTITUIR EL DESECANTE

Cuando abra el compartimento de desecante, impida la entrada de líquidos o gases inflamables en su interior. Para obtener instrucciones sobre el cambio de desecante, consulte <u>"Sustituir el desecante"</u>.

PELIGRO



Evite el peligro de explosión.

Antes de abrir el compartimento del desecante, apague el instrumento, desenchufe el cable de alimentación y extraiga las muestras del sistema. La entrada de líquidos o gases en el compartimento del desecante puede causar explosiones. Si esto ocurre, póngase en contacto con nosotros de inmediato y apague el instrumento hasta que se solucione el problema.

SUSTITUIR LA FUENTE IR

La fuente IR alcanza temperaturas extremadamente altas durante el funcionamiento. Si necesita sustituir la fuente IR, evite el peligro de quemaduras y explosiones.

PRECAUCIÓN



Evite el peligro de quemaduras.

La fuente IR alcanza temperaturas extremadamente altas durante el funcionamiento normal. Una vez apagado el espectrómetro, deje enfriar la fuente unos 10 minutos antes de trabajar con ella.

PELIGRO



Evite el peligro de explosión.

Antes de extraer la fuente del espectrómetro, apague el instrumento, desenchufe el cable de alimentación, desconecte las líneas de purga, y extraiga as muestras del sistema. La entrada de líquidos o gases en el compartimento de la fuente puede causar explosiones. Si esto ocurre, póngase en contacto con nosotros de inmediato y apague el instrumento hasta que se solucione el problema.

Resumen

El espectrómetro Nicolet Summit OA es un instrumento seguro y resistente, pero el usuario está expuesto a peligros durante su uso y mantenimiento. Durante el uso normal, tome precauciones cuando trabaje con muestras o disolventes peligrosos, y evite las partes del espectrómetro que están calientes o emiten calor. En las tareas de mantenimiento, tome precauciones para evitar lesiones o daños que pueden acontecer cuando se purga el instrumento, se manipulan componentes internos o se sustituye el desecante.

Desembalar y preparar el espectrómetro

Para desembalar y preparar adecuadamente el espectrómetro Nicolet Summit OA de Thermo Scientific necesita colocar el instrumento en el área de trabajo, conectar y encender el espectrómetro, y realizar una breve prueba de rendimiento.

Antes de comenzar

AVISO

Espere 24 horas antes de extraer el espectrómetro de la bolsa de plástico de transporte para evitar la formación de condensación, que puede dañar la óptica interna. La garantía no cubre daños causados por la apertura de la bolsa de plástico hermética antes de que el instrumento haya alcanzado la temperatura ambiente.

Recuerde: Repase la Información sobre seguridad y emplazamiento y asegúrese de la idoneidad y seguridad del área de trabajo escogido para el instrumento.

Desembalje y puesta en marcha

Los pasos para desembalar y preparar el espectrómetro varían en función de las opciones adquiridas con él.

1. Desembale la caja de transporte.

Extraiga el espectrómetro de la caja de transporte y la bolsa, y el resto de materiales de embalaje.

El espectrómetro lleva una tapa antipolvo extraíble sobre el cristal ATR. Utilice la tapa antipolvo cuando mueva o almacene el instrumento, y quítela cuando vaya a utilizar el instrumento.

- 2. Conecte el cable de alimentación al espectrómetro y a una fuente de corriente alterna debidamente conectada a tierra.
- 3. Conecte el monitor, el teclado y el ratón.
 - a. Conecte el monitor a la conexión Mini DisplayPort del espectrómetro.
 - b. Conecte el teclado y el ratón a uno de los puertos USB del espectrómetro.

Nota Se sabe que algunos dispositivos USB y Mini DisplayPort interfieren en la señal de dispositivos USB inalámbricos, como los teclados y ratones sin cables. Si observa problemas de funcionamiento con el teclado y el ratón inalámbricos, como retardos o pérdida de entradas, utilice el cable alargador USB provisto para colocar el receptor USB inalámbrico más alejado de los puertos USB.

4. Pulse el botón de alimentación del instrumento para encender el espectrómetro. La barra luminosa del espectrómetro muestra un ciclo de luz verde mientras el instrumento se calienta y una barra sólida de verde cuando el instrumento está listo para utilizarse.

Al encender o apagar el instrumento, puede oírse el sonido del movimiento de la referencia interna de poliestireno. Es normal.



Cuando se enciende el espectrómetro por primera vez, se le pide que introduzca una contraseña nueva. Le recomendamos que no introduzca ninguna contraseña en ese momento. Continúe sin introducir una contraseña y, una vez que haya terminado de configurarlo, gestione las cuentas y contraseñas del equipo de acuerdo con las pautas de su centro.

Nota Si la pantalla está en blanco al encender el sistema, compruebe si llega corriente al monitor, la conexión DisplayPort y la entrada del monitor. Consulte los detalles sobre el cambio de los ajustes de entrada en la documentación que acompaña al monitor.

5. Abra el software OMNIC Paradigm.

El software puede tardar varios minutos en iniciarse y conectarse al instrumento. El estado de conexión se muestra en el software en la esquina superior derecha de la pantalla. Cuando el estado indica que ya está conectado, continúe con el paso siguiente.



Nota Si abre cualquier menú del software antes de que el instrumento esté conectado, las opciones de alineación y calibración estarán desactivadas. Para activarlas, una vez que el instrumento esté conectado, haga clic en Panel de control, en la barra de herramientas.

- 6. Verifique el rendimiento del espectrómetro.
 - a. Alinee la óptica del espectrómetro.
 - Desde la interfaz de escritorio, vaya a Adquirir datos > Diagnósticos > Alinear espectrómetro.
 - ii. Seleccione **Iniciar** para comenzar la alineación. La pantalla muestra "Alineación completa" cuando termina de alinearse el instrumento.
 - b. Calibre el láser.
 - i. Vaya a Adquirir datos > Diagnósticos > Calibración de láser.
 - ii. Seleccione Iniciar para comenzar la calibración.
 - iii. La calibración del láser puede tardar varios minutos en realizarse. Una vez realizada la calibración, la pantalla muestra "Medición terminada" e indica cualquier cambio en la frecuencia del láser.

El instrumento necesita 12 horas para alcanzar el equilibrio térmico. Para unos resultados óptimos, repita la alineación y calibración después de 12 horas de funcionamiento.

c. Ejecute el diagrama de flujo Nicolet Summit OA - PV Test.

El diagrama de flujo de verificación del rendimiento (PV) emplea material de a referencia para realizar una serie de pruebas estándar y comprobar el funcionamiento del instrumento y la exactitud de los datos.

- i. Desde el panel·de·control de la interfaz de escritorio, desplácese hasta el panel de Diagramas de flujo.
- ii. En el panel de diagramas de flujo, haga clic con el botón derecho en el diagrama de flujo Nicolet Summit OA - PV Test y seleccione Ejecutar.

Primeros pasos con el espectrómetro 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. Siga las indicaciones de la pantalla para completar el diagrama de flujo.
- iv. Cuando se muestre la tabla de resultados de la prueba de verificación, seleccione Finalizar para completar el diagrama de flujo y volver al panel de control, o seleccione Measure Squalane Spectrum (Medir espectro de escualeno) para pasar a la sección de escualeno opcional del diagrama de flujo. La medición de un espectro de escualeno proporciona una prueba de referencia adicional para verificar el rendimiento del sistema.



Pasos siguientes

¡Enhorabuena! El instrumento ya está listo para utilizarse. No obstante, antes de comenzar la adquisición de datos, realice una copia de seguridad del sistema y cree una unidad de recuperación para proteger su sistema de problemas potenciales en el futuro. Windows 10 incluye varias herramientas internas para realizar copias de seguridad y recuperar datos.

Para obtener instrucciones sobre la creación de una unidad de recuperación e imagen del sistema, y sobre la restauración del sistema, consulte <u>"Copia de seguridad y recuperación de datos"</u>.

Si necesita más información sobre la identificación y el análisis de muestras con el espectrómetro, consulte las guías y los tutoriales de OMNIC Paradigm en thermofisher.com/ftir-help.

Para pasar a la interfaz de operador, vaya a Ver / Mostrar > Operador.

Encontrará instrucciones detalladas sobre la creación y ejecución de soluciones personalizadas con Package Editor y la interfaz Operator en "Custom Solutions", en la Guía del Usuario del software OMNIC Paradigm.

Copia de seguridad y recuperación de datos

Antes de comenzar la adquisición de datos, debe tomar algunas medidas de precaución para proteger su sistema de posibles problemas en el futuro. Windows 10 incluye varias herramientas internas para realizar copias de seguridad y recuperar datos.

¡IMPORTANTE!Para garantizar la recuperación de los datos en caso de que el disco duro o el sistema operativo resulten dañados, disponga de soportes de recuperación, cree una imagen del sistema y establezca procedimientos para realizar copias de seguridad de sus datos.

Crear una unidad de recuperación

Cree una unidad de recuperación para poder acceder a las opciones avanzadas de inicio de Windows 10 (Como Recuperación de imagen del sistema, Restauración del sistema y Configuración de inicio), incluso si su instalación de Windows resulta dañada.

Material necesario

• Unidad flash con formato FAT32 y un mínimo de 8 GB de espacio de almacenamiento

AVISO

En este proceso se eliminará todo el contenido de la unidad. Antes de continuar, asegúrese de haber realizado una copia de seguridad de todos los archivos que desee conservar.

Para crear una unidad de recuperación

- 1. Inserte la unidad flash en uno de los puertos USB del lateral del espectrómetro.
- 2. En el cuadro de búsqueda de la barra de tareas de Windows 10, busque "Crear una unidad de recuperación" y seleccione **Crear una unidad de recuperación**.



- 3. Seleccione Sí cuando se le pregunte si permite la realización de cambios en el equipo.
- 4. Anule la selección de "Realizar una copia de seguridad de los archivos de sistema en la unidad de recuperación" y haga clic en **Siguiente**.
- 5. Seleccione su unidad en la lista de unidades disponibles y haga clic en Siguiente.



- 6. Haga clic en **Crear**. Cuando haya terminado el proceso, haga clic en **Finalizar** para cerrar la ventana.
- 7. Retire con seguridad la unidad flash y guárdela en un lugar seguro. Una buena práctica es etiquetar la unidad como Unidad de Recuperación de Summit.

Para obtener más información sobre la creación de medios de recuperación, consulte la ayuda en línea de Microsoft.

Crear una imagen del sistema

Una imagen del sistema es una instantánea de todo el sistema, que incluye los datos, archivos, las aplicaciones instaladas y otros ajustes de configuración. En caso de fallo total del sistema operativo o del disco duro, puede utilizar la imagen del sistema para restaurar su equipo en el mismo estado en que estaba cuando se creó la imagen. Aunque es decisión suya cuándo actualizar la imagen del sistema, es recomendable conservar una imagen de la instalación limpia, tal como estuviera al configurar por primera vez el instrumento.

Este proceso es largo, aunque depende de la cantidad de datos que decida copiar en la imagen del sistema. Si se copia una gran cantidad datos, el proceso puede tardar más de una hora.

Material necesario

SSD con formato NTFS

Para crear una imagen del sistema

- 1. Conecte el SSD a uno de los puertos USB del espectrómetro.
- En Windows 10, abra el Panel de control y seleccione Copias de seguridad y restauración (Windows 7).
- 3. En el panel izquierdo, seleccione Crear una imagen de sistema.



4. Seleccione En un disco duro y elija su SSD en la lista. Haga clic en Siguiente.

| 🏥 Create a sys | tem image | | | |
|---|---|--|--|---------------------------------------|
| Where do y | ou want to save th | e backup? | | |
| A system image drives. A systen stops working; | is a copy of the drives i image can be used to however, you can't cho | required for Window restore your comput ose individual items | s to run. It can also ir er if your hard drive o to restore. | nclude additional or computer ever |
| On a hard di | sk | | | |
| - Backup | (E:) 465.60 GB free | | \sim | |
| | | | | |
| On one or m | ore DVDs | | _ | |
| On one or m | ore DVDs ve (D:) | | v | |
| ○ On one or m | ore DVDs ve (D:) k location | _ | × | |
| On one or m DVD Dri On a networ | ore DVDs ve (D:) k location | | Select | 1 |
| On one or m | ore DVDs ve (D:) k location | | ∽ Select | 1 |
| On one or m DVD Dri On a networ | ore DVDs ve (D:) k location | | ⊻ Select | I |

- 5. Confirme los ajustes de copia de seguridad y haga clic en Iniciar una copia de seguridad.
- 6. Responda No a la pregunta "¿Desea crear un disco de reparación del sistema?".
- 7. Haga clic en **Cerrar** y extraiga el SSD de forma segura del instrumento Summit. Etiquete la unidad como Recuperación de Summit y guárdela en un lugar seguro.

Restaurar el sistema

En caso de fallo del sistema operativo, utilice su unidad de recuperación e imagen del sistema para restaurar el sistema.

Material necesario

- Unidad USB de recuperación
- SSD con la imagen de su sistema
- Para restaurar el sistema con la unidad de recuperación y la imagen del sistema
- 1. Apague el espectrómetro Summit.
- 2. Conecte un teclado externo a uno de los puertos USB del espectrómetro Summit.
- 3. Conecte la unidad de recuperación a otro de los puertos USB.
- 4. Encienda el sistema.
- 5. Cuando aparezca la pantalla de presentación de Thermo Scientific, toque "**Esc**" en el teclado varias veces para abrir la BIOS del sistema.

Si aparece el cuadro de diálogo de "¿Desea salir sin guardar?" use las teclas de flecha para seleccionar "No".

- 6. Con las teclas de flecha, navegue hasta la pestaña Guardar y salir.
- 7. Navegue hasta la sección de anulación de arranque (Boot Override) y seleccione su unidad flash de recuperación. Pulse **Intro** para arrancar en la imagen de recuperación.



- 8. Cuando aparezca la pantalla "Elegir la distribución del teclado, extraiga del puerto USB la unidad flash de recuperación e inserte la SSD de imagen del sistema.
- 9. Seleccione la distribución de teclado que prefiera.
- 10. En la pantalla Elegir una opción, seleccione Solucionar problemas.

Primeros pasos con el espectrómetro Nicolet Summit OA

| Choo | ose an optio | n |
|---------------|---|---|
| \rightarrow | Continue Exit and continue to Windows 10 | |
| ľ | Troubleshoot Reset your PC or see advanced options | |
| Ċ | Turn off your PC | |
| | | |

- 11. Seleccione Opciones avanzadas.
- 12. En Opciones avanzadas, seleccione Recuperación de imagen del sistema.

| | System Restore Use a restore point recorded on your PC to restore Windows | < 0> | Startup Repair Fix problems that keep Windows from loading |
|----------|--|----------------|--|
| ₽ | Go back to the previous version | C:\ | Command Prompt Use the Command Prompt for advanced troubleshooting |
| _ | System Image Recovery Recover Windows using a specific system image the | ₿ | Startup Settings Change Windows startup behavior |

- 13. Seleccione Windows 10 como sistema operativo de destino.
- 14. En la ventana "Seleccionar una copia de seguridad de imagen del sistema", utilice la última imagen del sistema disponible o seleccione otra imagen del sistema. Haga clic en **Siguiente**.

| 🍋 Re-image your computer | | | × |
|---------------------------------|--|--|---|
| | Select a syst | em image backup | |
| | This computer will Everything on this information in the s | pe restored using the system image. computer will be replaced with the system image. | |
| | Troubleshooting in http://go.microsof | formation for BMR: t.com/fwlink/p/?LinkId=225039 | |
| | ◉ <u>U</u> se the latest a | vailable system image(recommended) | |
| | Location: | Backup (D:) | |
| | Date and time: | 9/3/2019 9:22:59 AM (GMT-8:00) | |
| | Computer: | DESKTOP-I6SNAFA | |
| ○ <u>S</u> elect a system image | | | |
| < Back Next > Cancel | | | |
| | | | _ |

15. Haga clic en **Siguiente** para dejar los valores predeterminados de las demás opciones de recuperación y haga clic en **Finalizar** para restaurar la imagen del sistema. Haga clic en **Yes** para confirmar.

La restauración tarda unos 15 minutos y, a continuación, el sistema se reinicia automáticamente.

Una vez recuperado el sistema, extraiga con seguridad la SSD del espectrómetro y guarde la unidad de recuperación y la SSD de imagen del sistema en un lugar seguro.

Para obtener más información sobre la restauración de la imagen del sistema, consulte la ayuda en línea de Microsoft.

Realizar copia de seguridad de los datos del sistema

Además de crear medios de recuperación y una imagen del sistema, es necesario que realice copias de seguridad de los datos de forma regular. Windows 10 incluye varias opciones para realizar copias de seguridad de los archivos, incluida la herramienta Historial de Archivos. El personal informático (TI) de su organización debe establecer los métodos para realizar las copias de seguridad más idóneos para las necesidades de su organización.

Instalar un adaptador inalámbrico USB

Puede conectar el espectrómetro Nicolet Summit OA a una red inalámbrica utilizando un adaptador inalámbrico USB opcional. Si se adquiere el adaptador con el espectrómetro, se suministra preinstalado.

Para instalar el adaptador inalámbrico USB, introdúzcalo en un puerto USB. Se desea realizar una instalación permanente, instale el adaptador en un puerto interior, bajo la cubierta del instrumento.

Para instalar un adaptador inalámbrico USB

1. Introduzca el adaptador en cualquier puerto USB disponible.



2. El sistema reconocerá el adaptador de forma automática. Para obtener información para conectarse a una red a través del adaptador, consulte la asistencia de Microsoft.

Para instalarlo en un puerto USB bajo la cubierta del instrumento

La instalación requiere quitar la placa de cristal y la cubierta del instrumento.

- 1. Siga las instrucciones de <u>"Sustituir la cubierta"</u> para quitar la placa de cristal y la cubierta del instrumento.
- 2. Cuando haya retirado la cubierta, introduzca el adaptador en uno de los puertos USB disponibles.



3. Vuelva a colocar la cubierta del instrumento y la placa de cristal.

Ahora el instrumento está listo para conectarse a una red inalámbrica.

Funcionamiento de la barra luminosa del espectrómetro

El espectrómetro Nicolet Summit OA de Thermo Scientific incluye una práctica barra luminosa que ofrece una rápida indicación visual de la calidad del producto y del estado del instrumento.

En las tablas siguientes se describen todas las señales de la barra luminosa.

CALENTAMIENTO EN CURSO

| Señal | Estado del sistema | Descripción |
|--------------|--|---|
| Vaivén verde | Sistema encendido y calentándose | La luz verde se desplaza de izquierda a derecha repetidamente. El tiempo aproximado de calentamiento es de 2 minutos. |

ADQUISICIÓN TRADICIONAL

| Señal | Estado del sistema | Descripción |
|--------------|--------------------------|---|
| Verde sólido | Preparado para el uso | Toda la barra luminosa se muestra en verde sólido. |
| Cometa verde | Adquiriendo datos | Sección corta de luz verde que oscila adelante y atrás. |

ADQUISICIÓN SMART

| Señal | Estado del sistema | Descripción |
|----------------|---|---|
| Pulsación azul | Adquiriendo fondo Smart | Toda la barra luminosa pulsa en azul. |
| Cometa azul | Adquiriendo datos de muestra (activado por usuario) | Sección corta de luz azul que oscila adelante y atrás. |

RESULTADOS DEL ANÁLISIS

| Señal | Estado del sistema | Descripción |
|----------------------|--|--|
| % relleno de verde | Valor de coincidencia o resultado de CCalidad por encima del umbral | El porcentaje de verde refleja el valor de coincidencia. Por ejemplo, un 90 % de relleno indica un valor de coincidencia de 90. |
| % relleno de naranja | Valor de coincidencia o resultado de CCalidad por debajo del umbral | El porcentaje de naranja refleja el valor de coincidencia. Por ejemplo, un 30 % de relleno indica un valor de coincidencia de 30. |

DIAGNÓSTICOS

| Señal | Estado del sistema | Descripción |
|-------------------|-----------------------|--|
| Destellos de rojo | Error del sistema | Toda la barra luminosa emite destellos rojos. Consulte la descripción del error en Estado del sistema , en el software OMNIC Paradigm. |

Ver los datos con la aplicación OMNIC Anywhere

OMNIC Anywhere es una aplicación basada en la nube que permite ver, examinar y compartir datos desde un PC, un ordenador Apple o un dispositivo Android o iOS.



Figura 2-1: OMNIC Anywhere en el navegador web

Con un espectrómetro Nicolet Summit OA y una cuenta Connect gratuita, puede medir muestras en el aula o el laboratorio, cargar los datos en su cuenta Connect y ver, examinar o compartir los datos con otro dispositivo en su residencia o área de trabajo.

Para obtener instrucciones sobre la creación de una cuenta Connect y ver sus datos con OMNIC Anywhere, visite "<u>Cloud-enabled FTIR Spectroscopy</u>" (Espectroscopia infrarroja por transformada de Fourier (FTIR) en la nube). [Página en blanco]

Capítulo 3

Opciones y accesorios

El Nicolet Summit OA se puede utilizar con el kit de purga opcional.

| nstalar y mantener un kit de purga34 |
|--------------------------------------|
|--------------------------------------|

Instalar y mantener un kit de purga

La purga del espectrómetro con aire seco o nitrógeno protege los componentes internos de la humedad y otros contaminantes ambientales. Para instalar el kit de purga necesita montar las válvulas y los reguladores, sustituir el cartucho de desecante, conectar el gas de purga y establecer la presión y el caudal.

ADVERTENCIA



Evite el riesgo de explosión.

No utilice nunca un gas tóxico, inflamable o combustible para purgar el instrumento. El gas de purga debe estar libre de aceite y otros materiales reactivos. El calor producido por la fuente o por la absorción de láser puede encender los gases inflamables o los materiales reactivos que se encuentran en el gas de purga. Use solo nitrógeno o aire seco para purgar el instrumento.

AVISO

Se recomienda mantener el instrumento sellado y desecado o purgado en todo momento. La garantía no cubre daños por la falta de desecación o purga adecuadas del instrumento. Si tiene alguna duda a este respecto, póngase en contacto con nosotros.

CONFIGURACIÓN NECESARIA

Antes de instalar el kit de purga, necesita contar con una fuente de aire seco o nitrógeno que cumpla las características técnicas especificadas por Thermo Scientific para el espectrómetro Nicolet[™] Summit. Consulte los detalles en la Guía de seguridad y emplazamiento del instrumento.

Nota Para obtener los mejores resultados, seque el gas de purga hasta un punto de rocío de -70 °C (-94 °F) o menos.

MATERIAL NECESARIO

Además del kit de purga, necesita el material y las herramientas siguientes:

- Llave fija abierta de 3/4 pulgada
- Llave fija abierta de 11/16 pulgada
- Destornillador con punta Phillips n.º 2
- Llave hexagonal de 2,381 mm (3/32 pulg.)

Opciones y accesorios

- Llave hexagonal de 2,778 mm (7/64 pulg.)
- Cinta de sellado de roscas ("cinta de fontanero" o "cinta de teflón")

Para instalar un kit de purga

- 1. Conecte el conjunto de fontanería a la fuente de gas de purga.
 - a. Instale la válvula reguladora y un racor macho de 1/4 pulgada o uno hebra de 3/8 pulgada en la fuente de gas de purga. (Elija una válvula y racores que sean apropiados para la fuente de gas de purga).



b. Si utiliza un racor macho de 1/4 pulgada, continúe en el paso siguiente.

Si utiliza un racor hembra de 3/8 pulgada en la fuente de gas de purga, debe instalar la boquilla reductora de 3/8 a 1/4 pulgada incluida en el kit de purga. Aplique cinta de teflón a la boquilla reductora antes de instalarla, y utilice una llave fija de boca abierta de 11/16 pulgada para apretar la conexión.



Opciones y accesorios

c. Aplique cinta de teflón a la boquilla reductora o al racor macho de 1/4 pulgada e instale el acoplamiento de presión. Utilice una llave fija de boca abierta de 3/4 pulgada para apretar la conexión.



- d. Encaje con firmeza la entrada macho del conjunto de fontanería de la pared en el racor de desenganche rápido.
- 2. Apague el espectrómetro.
 - a. Para apagar el espectrómetro, pulse el botón de encendido/apagado.
 - b. Desenchufe el cable de alimentación.
- 3. "Quitar la placa de cristal"
- 4. Retire la cubierta del instrumento.
 - a. Incline con cuidado el instrumento hacia atrás hasta que quede apoyado en las aletas del disipador de calor.
 - b. Con la llave hexagonal 7/64, afloje los cuatro tornillos que fijan la cubierta de la base del instrumento.



- c. Con cuidado, vuelva a colocar el instrumento sobre las patas. Al enderezar el instrumento, se caerán los cuatro tornillos. Tenga cuidado de no perderlos.
- d. Levante la cubierta hacia arriba en vertical desde la base.



5. Extraiga el cartucho de desecante.

Opciones y accesorios

a. Utilice el destornillador con punta Phillips n.º 2 para aflojar los dos tornillos cautivos del cartucho de desecante y levante el cartucho recto hacia arriba para extraerlo del instrumento.



b. Extraiga y deseche la junta tórica grande.



Nota Tenga cuidado de que no caiga la junta tórica dentro del compartimento de desecante.

- 6. Abra el kit de purga y el paquete de desecante sellado.
- 7. Extraiga la junta tórica del paquete y colóquela en la placa de asiento del instrumento (ver imagen anterior). Presione (mejor que enrollarla) la nueva junta tórica en su lugar para asegurarse de que queda bien encajada en la ranura.

Opciones y accesorios

8. Extraiga el cartucho de desecante con racores de purga del paquete, inserte el cartucho en el compartimento de desecante, compruebe que queda bien asentado en la junta tórica y fíjelo mediante los dos tornillos.

Fíjese en la orientación correcta del cartucho. Cuando el cartucho está instalado, se puede leer su etiqueta desde delante del instrumento.



- 9. Dirija la línea de purga por la apertura de entrada de purga, en la parte trasera del instrumento.
 - a. Baje con cuidado la cubierta sobre la base mientras coloca el tubo de purga en la ranura de la cubierta.



- b. Baje la cubierta completamente. Al bajar la cubierta, el tubo de purga queda atrapado en su ranura entre la cubierta y el disipador de calor.
- c. Aparte el tubo de purga a un lado e incline el instrumento hacia atrás, de modo que quede apoyado en las aletas del disipador de calor.
- d. Inserte los cuatro tornillos que retiró previamente. Apriete los tornillos.
- e. Vuelva a colocar el instrumento sobre las patas.
- f. Introduzca el exceso de tubo en el instrumento, o tire de él, según proceda.



10. <u>"Volver a colocar la placa de cristal"</u>.

- 11. Conecte el acoplador de flujo del conjunto de fontanería de la pared (marcado con "To spectrometer" [Al espectrómetro]) al tubo de purga.
- 12. Ajuste los controles del gas de purga.
 - a. Abra la válvula de regulación principal y ajuste los controles del gas de purga como se indican a continuación (consulte Establecer controles del gas de purga para más detalles).

| Hardware | Presión (psig) | Caudal (scfh) |
|-------------------------|----------------|---------------|
| Espectrómetro Summit | 5 | 1 |

- b. Enchufe el cable de alimentación al instrumento y enciéndalo.
- c. Espere 30-60 minutos para que el instrumento se purgue completamente.

Establecer controles del gas de purga

El ajuste adecuado de los controles del gas de purga garantiza la protección antihumedad del espectrómetro sin introducir vibraciones.

Para unos resultados óptimos, el gas de purga se debe secar hasta un punto de rocío de - 70ºC (- 94ºF) o inferior.



- 1. Abra la válvula principal para que comience a circular gas de purga por el regulador.
- 2. Ajuste el regulador de presión del espectrómetro hasta que el manómetro indique una presión de 5 psig (34 kPa).
- 3. Ajuste el caudalímetro del espectrómetro a 1 scfh (0,47 l/min).

AVISO

Caudales superiores a los recomendados pueden originar vibraciones y afectar a la calidad de los datos.

Comprobar y cambiar el filtro del gas de purga

Sustituya el filtro de purga cuando esté amarillo o descolorido, o si muestra restos de suciedad o partículas extrañas.

4.

AVISO

Se recomienda mantener el instrumento sellado y desecado o purgado en todo momento. Los daños causados al equipo por no mantener el instrumento sellado y desecado, o sellado y purgado, no están cubiertos por la garantía. Si tiene alguna duda a este respecto, póngase en contacto con nosotros.

Figura 3-1: Filtro de purga dentro de la cazoleta de plástico situada bajo el manómetro



Nota Para pedir componentes, póngase en contacto con Soporte Técnico.

Para cambiar el filtro del gas de purga

- 1. Cierre el gas de purga en la válvula principal. No baje el caudalímetro ni el regulador de presión.
- 2. Retire la cazoleta de plástico que contiene el filtro y extraiga el filtro. (Puede desenroscarlos con la mano).



Opciones y accesorios

- 3. Introduzca el nuevo filtro y vuelva a instalar la cazoleta.
- 4. Abra la válvula principal para que pase el gas de purga al instrumento y compruebe la corrección del caudal de gas de purga.

[Página en blanco]

Capítulo 4

Mantenimiento

El espectrómetro FTIR Nicolet Summit OA requiere muy pocos cuidados y ningún mantenimiento diario. No obstante, necesita de los cuidados básicos para mantener su máximo rendimiento.

| Programa de mantenimiento del espectrómetro Summit | 46 |
|--|----|
| Alinear la óptica del espectrómetro | |
| Calibrar la frecuencia del láser | |
| Limpiar el instrumento y el cristal | |
| Quitar y sustituir la placa de cristal | 52 |
| Sustituir la cubierta | 54 |
| Sustituir el desecante | |
| Sustituir la fuente | 64 |

Programa de mantenimiento del espectrómetro Summit

El espectrómetro FTIR Nicolet Summit OA de Thermo Scientific requiere muy pocos cuidados y ningún mantenimiento diario. No obstante, necesita de los cuidados básicos para mantener su máximo rendimiento. Siga las pautas de este artículo. Para un rendimiento óptimo, deje el espectrómetro encendido.

AVISO

La electricidad estática puede dañar irreversiblemente los componentes esenciales del espectrómetro. Para evitar tales daños, siga estas recomendaciones:

- Antes de desconectar la fuente de alimentación, toque la base metálica del espectrómetro para descargar cualquier electricidad estática que usted pueda tener acumulada.
- Deje las piezas de sustitución en su envase protector hasta que esté preparado para instalarlas en el instrumento.

Mantenimiento semanal

VERIFICAR EL RENDIMIENTO DEL ESPECTRÓMETRO

El software OMNIC Paradigm incluye un diagrama de flujo de verificación del rendimiento (PV) para el espectrómetro. El diagrama de flujo PV ejecuta una serie de pruebas estándar para comprobar el funcionamiento del instrumento y garantizar la exactitud de los datos.

Recomendamos ejecutar el diagrama de flujo PV al menos una vez a la semana.

LIMPIAR EL ESPECTRÓMETRO

El espectrómetro se debe limpiar solo según las recomendaciones descritas. Consulte <u>"Limpiar el</u> espectrómetro".

Mantenimiento mensual

COMPROBAR EL INDICADOR DE HUMEDAD

Los componentes ópticos del espectrómetro se pueden dañar fácilmente por el exceso de humedad en el aire. El espectrómetro está sellado y sus componentes está protegidos por dos recipientes de

desecante que absorben la humedad. Controle la humedad del instrumento en el software OMNIC Paradigm.

AVISO

Se recomienda mantener el espectrómetro sellado y desecado, o sellado y purgado con aire seco no nitrógeno, en todo momento. Los daños causados al equipo por no mantener el instrumento sellado y desecado, o sellado y purgado, no están cubiertos por la garantía. Si tiene alguna duda a este respecto, póngase en contacto con nosotros.

COMPROBAR EL FILTRO DEL GAS DE PURGA

Si el espectrómetro se purga con nitrógeno o aire seco, compruebe el filtro de purga al menos una vez al mes.

Sustituya el filtro purga cuando esté amarillo o descolorido, o si muestra restos de suciedad o partículas extrañas. Para obtener más información, consulte "Comprobar y cambiar el filtro del gas de purga".

Figura 4-1: Filtro de purga dentro de la cazoleta de plástico situada bajo el manómetro



Alinear la óptica del espectrómetro

Alinee la óptica del espectrómetro si el diagrama de flujo Nicolet Summit OA - PV Test indica que no ha superado la prueba. Alinear el espectrómetro optimiza la energía que llega al detector y maximiza su señal. Si no se supera la prueba de rendimiento, alinee la óptica del espectrómetro, calibre la frecuencia del láser y vuelva a realizar la prueba. Si continúa sin superarse la prueba, debe solicitar ayuda al representante local de Soporte Técnico.

- Para alinear la óptica del espectrómetro (interfaz de pantalla táctil)
- 1. En la pantalla de inicio, seleccione el icono Diagnósticos [200] para abrir la vista de diagnósticos.
- 2. Abra la pestaña Alinear y seleccione Alinear.

El sistema alinea automáticamente el espectrómetro e informa de su finalización.

- Para alinear la óptica del espectrómetro (interfaz de operador)
- 1. Desplácese hasta el área del Administrador y seleccione el icono Diagnósticos [
- 2. Abra la pestaña Alinear y seleccione Alinear.

El sistema alinea automáticamente el espectrómetro e informa de su finalización.

- Para alinear la óptica del espectrómetro (interfaz de escritorio)
- 1. Seleccione Adquirir datos > Diagnósticos > Alinear espectrómetro.
- 2. En el cuadro de diálogo Alinear, haga clic en Inicio.

El sistema alinea automáticamente el espectrómetro e informa de su finalización.

Calibrar la frecuencia del láser

Calibre la frecuencia del láser si el diagrama de flujo Nicolet Summit OA - PV Test indica que no ha superado la prueba. Si no se supera una prueba, es preciso alinear primero el espectrómetro, calibrar después la frecuencia del láser y, a continuación, volver a realizar la prueba. Si continúa sin superarse la prueba, debe solicitar ayuda al representante local de Soporte Técnico.

Para calibrar el láser (interfaz de pantalla táctil)

- 1. En la pantalla de inicio, seleccione el icono Diagnósticos [
- 2. Abra la pestaña Calibración de láser y seleccione Calibrar para comenzar.

Cuando finaliza la calibración del láser, un mensaje indica si la frecuencia ha cambiado e informa de la nueva frecuencia.

Para calibrar el láser (interfaz de operador)

- 1. Desplácese hasta el área del Administrador y seleccione el icono Diagnósticos [
- 2. Abra la pestaña Calibración de láser y seleccione Calibrar para comenzar.

Cuando finaliza la calibración del láser, un mensaje indica si la frecuencia ha cambiado e informa de la nueva frecuencia.

Para calibrar el láser (interfaz de escritorio)

- 1. Seleccione Adquirir datos > Diagnósticos > Calibración de láser.
- 2. En el cuadro de diálogo Calibración de láser, haga clic en Inicio para calibrar el láser.

Cuando el sistema termina de calibrar el láser, muestra un mensaje indicando si la frecuencia ha cambiado e informa de la nueva frecuencia.

Limpiar el instrumento y el cristal

Antes de limpiar el espectrómetro Thermo Scientific Nicolet Summit OA, apáguelo y desconecte el cable de alimentación.

Limpiar el espectrómetro

Limpie con cuidado el exterior del espectrómetro con un paño suave limpio y un disolvente suave. Rocíe el paño y limpie exclusivamente la superficie pintada exterior. Tenga cuidado y evite que el disolvente llegue a los componentes electrónicos de la parte trasera del instrumento.

Podría depositarse polvo en el recinto de componentes electrónicos, en la parte trasera del instrumento, e interferir en la disipación de calor, lo que reduciría la vida útil de dichos componentes.

Para eliminar el polvo de la parte trasera del instrumento, sóplelo con aire comprimido. No utilice líquido alguno para quitar el polvo de la parte trasera del espectrómetro.

PRECAUCIÓN



Evite el peligro de descargas eléctricas. Para evitar el peligro de descargas, no permita que entre líquido en la fuente de alimentación ni en la parte trasera del instrumento.

AVISO

No use detergentes fuertes, disolventes ni substancias químicas o abrasivas porque pueden dañar el acabado del instrumento.

Limpiar la placa del cristal ATR

Para limpiar el cristal ATR, utilice un paño o pañuelo limpio, seco y no abrasivo.

ADVERTENCIA



Evite la inhalación e ingestión de sustancias tóxicas.

El seleniuro de zinc (ZnSe) es tóxico. Consulte las medidas de control de manipulación y exposición en la Ficha de datos de seguridad de ZnSe del fabricante, en www.specac.com.

Lleve guantes protectores cuando limpie o sustituya la placa de cristal.

- 1. Seque el cristal con un paño o tejido. Deseche el paño usado.
- 2. Con una pipeta de transferencia, aplique algunas gotas de heptano al cristal y sus alrededores.
- 3. Pase un paño limpio nuevo a lo largo del cristal para retirar cualquier rastro de líquido. Limpie también el área en torno al cristal.



4. Repita la operación, sin es necesario, acabando siempre con un paño limpio y seco.

Quitar y sustituir la placa de cristal

La placa de cristal se debe sustituir cada 12-18 meses, según sea necesario.

ADVERTENCIA



Evite la inhalación e ingestión de sustancias tóxicas.

El seleniuro de zinc (ZnSe) es tóxico. Consulte las medidas de control de manipulación y exposición en la Ficha de datos de seguridad de ZnSe del fabricante, en www.specac.com.

Lleve guantes protectores cuando limpie o sustituya la placa de cristal.

AVISO

Tenga cuidado y evite que caiga cualquier cosa en el instrumento cuando haya quitado la cubierta.

MATERIAL NECESARIO

• Llave hexagonal de 2,381 mm (3/32 pulg.)

Quitar la placa de cristal

- 1. Retire la tapa antipolvo del cristal.
- 2. Con una llave hexagonal de 3/32 pulg. (2,381 mm), afloje totalmente los tres tornillos que sujetan la placa de cristal al instrumento.



- 3. Levante la placa de cristal hacia arriba en vertical desde el instrumento.
- 4. Utilice un paño limpio y seco para limpiar el área alrededor del reborde, debajo de la zona de asiento de la placa.

Volver a colocar la placa de cristal

- 1. Ajuste con precisión la nueva placa sobre el reborde.
- 2. Coloque las arandelas en los tornillos y suelte los tornillos en los tres orificios. Los tornillos deben colocarse dentro de los orificios.
- Mientras presiona la placa de cristal, apriete los dos tornillos de los lados de la placa y, a continuación, el tercer tornillo, más próximo a la parte trasera del instrumento. Apriete todos los tornillos a mano.

Si el tornillo queda suelto, desenrósquelo y vuelva a insertarlo para corregir su alineación.

Sustituir la cubierta

HERRAMIENTAS NECESARIAS

Antes de comenzar, compruebe que cuenta con lo siguiente:

- Llave hexagonal 7/64
- Llave hexagonal 3/32

Para sustituir la cubierta del espectrómetro

- 1. Apague el espectrómetro.
 - a. Para apagar el espectrómetro, pulse el botón de encendido/apagado.
 - b. Desenchufe el cable de alimentación.
 - c. Desactive la purga, si procede, y desconecte la línea de purga que llega al instrumento.
- 2. Retire la placa de cristal.
 - a. Retire la tapa antipolvo del cristal.
 - b. Afloje totalmente los tres tornillos que sujetan la placa de cristal al instrumento.



- c. Levante la placa de cristal hacia arriba en vertical desde el instrumento.
- d. Utilice un paño limpio y seco para limpiar el área alrededor del reborde, debajo de la zona de asiento de la placa.
- 3. Retire la cubierta del instrumento.
 - a. Incline con cuidado el instrumento hacia atrás hasta que quede apoyado en las aletas del disipador de calor.
 - b. Con la llave hexagonal 7/64, afloje los cuatro tornillos que fijan la cubierta de la base del instrumento.



c. Con cuidado, vuelva a colocar el instrumento sobre las patas. Al enderezar el instrumento, se caerán los cuatro tornillos. Tenga cuidado de no perderlos.

d. Levante la cubierta hacia arriba en vertical desde la base.



- 4. Vuelva a colocar la cubierta del instrumento.
 - a. Ajuste con precisión la cubierta en la base del instrumento.
 - b. Incline el instrumento hacia atrás apoyándolo en las aletas del disipador de calor.
 - c. Inserte los cuatro tornillos que retiró previamente. Apriete los tornillos.
 - d. Vuelva a colocar el instrumento sobre las patas.
- 5. Vuelva a colocar la placa de cristal.
 - a. Ajuste con precisión la placa sobre el reborde.
 - b. Coloque las arandelas en los tornillos y suelte los tornillos en los tres orificios. Los tornillos deben colocarse dentro de los orificios.
 - c. Mientras presiona la placa de cristal, apriete los dos tornillos de los lados de la placa y, a continuación, el tercer tornillo, más próximo a la parte trasera del instrumento. Apriete todos los tornillos a mano.

Si el tornillo queda suelto, desenrósquelo y vuelva a insertarlo para corregir su alineación.

6. Encienda el espectrómetro.

- a. Enchufe el cable de alimentación al instrumento y enciéndalo.
- b. Vuelva a conectar la línea de purga al instrumento y, si procede, active la purga.
- 7. Verifique el rendimiento del espectrómetro.
 - a. Alinee la óptica del espectrómetro.
 - Desde el panel de control del software OMNIC Paradigm, vaya a Adquirir datos
 > Diagnósticos > Alinear espectrómetro.
 - ii. Seleccione **Iniciar** para comenzar la alineación. La pantalla muestra "Alineación completa" cuando termina de alinearse el instrumento.
 - b. Calibre el láser.
 - i. Vaya a Adquirir datos > Diagnósticos > Calibración de láser.
 - ii. Seleccione Iniciar para comenzar la calibración.

La calibración del láser puede tardar varios minutos en realizarse. Una vez realizada la calibración, la pantalla muestra "Medición terminada" e indica cualquier cambio en la frecuencia del láser.

El instrumento necesita 12 horas para alcanzar el equilibrio térmico. Para unos resultados óptimos, repita la alineación y calibración después de 12 horas de funcionamiento.

- c. Ejecute el diagrama de flujo Nicolet Summit OA PV Test. El diagrama de flujo de verificación del rendimiento (PV) emplea material de referencia para realizar una serie de pruebas estándar y comprobar el funcionamiento del instrumento y la exactitud de los datos.
 - i. Vuelva al panel de control.
 - ii. Desplácese hasta el panel de Diagramas de flujo y seleccione Nicolet Summit OA -PV Test.
 - iii. Haga clic con el botón derecho del ratón y seleccione **Ejecutar** para comenzar el diagrama de flujo.

Sustituir el desecante

Controle la humedad del instrumento mediante el software OMNIC Paradigm. Sustituya el desecante cuando sea necesario o según las pautas del centro.

Para cambiar el desecante es preciso retirar la cubierta del instrumento, extraer el cartucho de desecante y sustituir la junta tórica y los recipientes de desecante.

HERRAMIENTAS NECESARIAS

Antes de comenzar, compruebe que cuenta con lo siguiente:

- Llave hexagonal 7/64
- Kit de sustitución de desecante
- Llave hexagonal 3/32
- Destornillador con punta Phillips n.º 2
- Guantes, dediles o paño de laboratorio (para manipular el indicador de humedad)

Evite el riesgo de explosión.

ADVERTENCIA



Antes de retire la cubierta del instrumento y abrir el compartimento del desecante, apague el instrumento, desenchufe el cable de alimentación y extraiga todas las muestras del sistema.

La entrada de líquidos o gases en el compartimento de desecante conlleva peligro de explosiones. Si esto ocurre, póngase en contacto con nosotros de inmediato y apague el instrumento hasta que se solucione el problema.

AVISO

Asegúrese de que no caiga nada dentro del instrumento mientras está retirada la cubierta del desecante.

ADVERTENCIA



Cuando sustituya el desecante, utilice solo las piezas de repuesto suministradas por nosotros.

Para cambiar el desecante

- 1. Apague el espectrómetro.
 - a. Para apagar el espectrómetro, pulse el botón de encendido/apagado.
 - b. Desenchufe el cable de alimentación.
 - c. Desactive la purga, si procede, y desconecte la línea de purga que llega al instrumento.

Nota Para mantener la humedad requerida, puede utilizar cualquiera de los cartuchos de desecante (con o sin conector de purga). Las siguientes ilustraciones muestran la versión de cartucho de desecante no purgado. Las instrucciones sobre el cambio de recipientes de desecante son las mismas para ambos cartuchos.

- 2. "Quitar la placa de cristal"
- 3. Retire la cubierta del instrumento.
 - a. Incline con cuidado el instrumento hacia atrás hasta que quede apoyado en las aletas del disipador de calor.
 - b. Con la llave hexagonal 7/64, afloje los cuatro tornillos que fijan la cubierta de la base del instrumento.



c. Con cuidado, vuelva a colocar el instrumento sobre las patas. Al enderezar el instrumento, se caerán los cuatro tornillos. Tenga cuidado de no perderlos.

d. Levante la cubierta hacia arriba en vertical desde la base.



- 4. Extraiga el cartucho de desecante.
 - a. Utilice el destornillador con punta Phillips n.º 2 para aflojar los dos tornillos cautivos del cartucho de desecante y levante el cartucho recto hacia arriba para extraerlo del instrumento.



b. Extraiga y deseche la junta tórica grande.



Nota Tenga cuidado de que no caiga la junta tórica dentro del compartimento de desecante.

5. Extraiga los recipientes de desecante y sustituya el indicador de humedad.

ADVERTENCIA



El contenido de los recipientes de desecante (tamiz molecular de aluminosilicato metálico hidratado) puede ser nocivo si se ingiere. Cuando deseche los recipientes de desecante saturado, asegúrese de eliminarlos de la forma adecuada.

a. Utilice la llave hexagonal de 1,27 mm para aflojar (unas dos vueltas) los ocho tornillos que sujetan los recipientes de desecante saturado al cartucho de desecante. Extraiga los recipientes.



b. Gire hacia abajo el cartucho de desecante y despegue del cristal el indicador de humedad antiguo. Deseche el indicador usado.

AVISO

Use siempre guantes, dediles o paño de laboratorio cuando manipule el indicador de humedad nuevo. La grasa o humedad de la piel puede decolorar el indicador.

- 6. Instale los nuevos recipientes de desecante.
 - a. Abra el paquete sellado que contiene el desecante nuevo (debe estar precintado para conservar bien el desecante) y extraiga en indicador de humedad nuevo.
 - b. Gire hacia abajo el cartucho de desecante y presione el indicador azul sobre el cristal para que los bordes planos del indicador queden al nivel de los bordes del cartucho. El texto debe quedar visible a través del cristal cuando vuelva a colocar el cartucho hacia arriba.

AVISO

Asegúrese de que el indicador quede bien adherido al cristal para que no se suelte del compartimento de desecante al volver a colocar el cartucho.

- c. Inserte los recipientes en el canal del cartucho de desecante y apriete los ocho tornillos (unas 2 vueltas) hasta que queden alineados en el canal del cartucho.
- 7. Introduzca el cartucho de desecante.
 - a. Extraiga la junta tórica nueva del paquete y colóquela en la placa de asiento del instrumento.

Presione (mejor que enrollarla) la nueva junta tórica en su lugar para asegurarse de que queda bien encajada en la ranura.

b. Introduzca el cartucho de desecante en el espectrómetro de forma que el texto del indicador de humedad pueda leerse estando delante del instrumento.

Asegúrese de que el cartucho quede bien asentado sobre la junta tórica y apriete los dos tornillos cautivos con el destornillador de punta Phillips n.º 2.

- 8. Vuelva a colocar la cubierta del instrumento.
 - a. Ajuste con precisión la cubierta en la base del instrumento.
 - b. Incline el instrumento hacia atrás apoyándolo en las aletas del disipador de calor.
 - c. Inserte los cuatro tornillos que retiró previamente. Apriete los tornillos.

- d. Vuelva a colocar el instrumento sobre las patas.
- e. "Volver a colocar la placa de cristal".
- 9. Encienda el espectrómetro.
 - a. Enchufe el cable de alimentación al instrumento y enciéndalo.
 - b. Vuelva a conectar la línea de purga al instrumento y, si procede, active la purga.
- 10. Verifique el rendimiento del espectrómetro.
 - a. Alinee la óptica del espectrómetro.
 - i. Desde el panel de control del software OMNIC Paradigm, vaya a Adquirir datos
 > Diagnósticos > Alinear espectrómetro.
 - ii. Seleccione **Iniciar** para comenzar la alineación. La pantalla muestra "Alineación completa" cuando termina de alinearse el instrumento.
 - b. Calibre el láser.
 - i. Vaya a Adquirir datos > Diagnósticos > Calibración de láser.
 - ii. Seleccione Iniciar para comenzar la calibración.

La calibración del láser puede tardar varios minutos en realizarse. Una vez realizada la calibración, la pantalla muestra "Medición terminada" e indica cualquier cambio en la frecuencia del láser.

El instrumento necesita 12 horas para alcanzar el equilibrio térmico. Para unos resultados óptimos, repita la alineación y calibración después de 12 horas de funcionamiento.

- c. Ejecute el diagrama de flujo Nicolet Summit OA PV Test. El diagrama de flujo de verificación del rendimiento (PV) emplea material de referencia para realizar una serie de pruebas estándar y comprobar el funcionamiento del instrumento y la exactitud de los datos.
 - i. Vuelva al panel de control.
 - ii. Desplácese hasta el panel de Diagramas de flujo y seleccione Nicolet Summit OA -PV Test.
 - iii. Haga clic con el botón derecho del ratón y seleccione **Ejecutar** para comenzar el diagrama de flujo.

Sustituir la fuente

La fuente IR es fácilmente accesible desde la base del espectrómetro y se puede cambiar sin necesidad de retirar la cubierta del espectrómetro.

HERRAMIENTAS Y MATERIAL NECESARIOS

- Destornillador con punta Phillips n.º 1
- Kit de sustitución de fuente

PRECAUCIÓN



Antes de sustituir la fuente, apague el instrumento. Cuando sustituya la fuente, utilice solo las piezas de repuesto suministradas por nosotros.

Para cambiar la fuente

- 1. Apague y desenchufe el espectrómetro.
 - a. Apague el instrumento y desenchufe el cable de alimentación.
 - b. Desconecte todos los cables del instrumento, incluidos el cable Ethernet y los dispositivos USB.
 - c. Si el espectrómetro está purgado, desconectéctele las líneas de purga. Consulte <u>"Instalar y</u> <u>mantener un kit de purga"</u> para obtener más información.

PRECAUCIÓN



d.

Evite el peligro de quemaduras.

La fuente alcanza temperaturas extremadamente altas durante el funcionamiento del instrumento. Tras apagar el instrumento, espere al menos 10 minutos antes de realizar el paso siguiente.

- 2. Extraiga la fuente.
 - a. Utilice el destornillador con punta Phillips n.º 1 para aflojar completamente los tres tornillos de sujeción de la fuente.



b. Gire y tire de la placa hacia fuera. Sujete el instrumento con una mano y tire de la placa hacia fuera con decisión para extraer la fuente del instrumento.



AVISO

En ningún momento debe tocar la resistencia de la fuente directamente con los dedos. El aceite de la piel u otros residuos acortan la vida útil de la resistencia. Utilice siempre guantes, dediles o paños de laboratorio limpios cuando manipule la resistencia de una fuente.

- 3. Inserte la fuente nueva.
 - a. Asegúrese de que los tres tornillos cautivos estén derechos y, a continuación, introduzca la fuente en el compartimento hasta que el conjunto de la fuente esté alineado con el instrumento.
 - b. Apriete los tornillos para fijar la fuente y encaje la placa en la su posición.
- 4. Vuelva a enchufar y a encender el espectrómetro.
 - a. Coloque de nuevo el instrumento en posición vertical lentamente y vuelva conectar los cables retirados con anterioridad.

b. Sustituya el desecante. Consulte las instrucciones en "Sustituir el desecante".

Debido a la exposición de la óptica interna a la atmósfera al retirar la fuente, siempre es preciso cambiar el desecante después de sustituir la fuente.

ADVERTENCIA



Evite riesgos de descarga eléctrica.

Para asegurar una buena conexión a tierra y evitar el peligro de descargas, no utilice tomas de corriente que estén conectadas a un conducto de tierra. La toma de tierra debe ser un cable que no lleve corriente y que esté conectado al punto de tierra en la caja principal de distribución.

- c. Enchufe el cable de alimentación al instrumento y pulse el botón de encendido/apagado para encender el espectrómetro.
- 5. Verifique el rendimiento del espectrómetro.
 - a. Alinee la óptica del espectrómetro.
 - Desde el panel de control del software OMNIC Paradigm, vaya a Adquirir datos
 > Diagnósticos > Alinear espectrómetro.
 - ii. Seleccione **Iniciar** para comenzar la alineación. La pantalla muestra "Alineación completa" cuando termina de alinearse el instrumento.
 - b. Calibre el láser.
 - i. Vaya a Adquirir datos > Diagnósticos > Calibración de láser.
 - ii. Seleccione Iniciar para comenzar la calibración.

La calibración del láser puede tardar varios minutos en realizarse. Una vez realizada la calibración, la pantalla muestra "Medición terminada" e indica cualquier cambio en la frecuencia del láser.

El instrumento necesita 12 horas para alcanzar el equilibrio térmico. Para unos resultados óptimos, repita la alineación y calibración después de 12 horas de funcionamiento.

c. Ejecute el diagrama de flujo Nicolet Summit OA - PV Test. El diagrama de flujo de verificación del rendimiento (PV) emplea material de referencia para realizar una serie de pruebas estándar y comprobar el funcionamiento del instrumento y la exactitud de los datos.

- i. Vuelva al panel de control.
- ii. Desplácese hasta el panel de Diagramas de flujo y seleccione Nicolet Summit OA -PV Test.
- iii. Haga clic con el botón derecho del ratón y seleccione **Ejecutar** para comenzar el diagrama de flujo.

[Página en blanco]