Nicolet Summit OA

FTIR Spektrometer



BEDIENUNGSANLEITUNG

269-346500 Ausgabe B April 2021

© 2020 Thermo Fisher Scientific Inc. Alle Rechte vorbehalten.

Wenden Sie sich für einen technischen Support an www.thermofisher.com.

Dieses Dokument liegt allen Produkten von Thermo Fisher Scientific Inc. beim Kauf bei und ist beim Betrieb des Produkts zu beachten. Dieses Dokument ist urheberrechtlich geschützt. Jede teilweise oder vollständige Kopie dieses Dokuments ist streng untersagt, sofern keine schriftliche Genehmigung von Thermo Fisher Scientific Inc. vorliegt.

Der Inhalt dieses Dokuments kann jederzeit ohne Ankündigung geändert werden. Sämtliche technische Informationen in diesem Dokument dienen lediglich zu Referenzzwecken. In diesem Dokument genannten Systemkonfigurationen und -spezifikationen ersetzen alle vorher gegebenen Informationen.

Thermo Fisher Scientific Inc. erhebt keinen Anspruch auf Vollständigkeit, Genauigkeit und Fehlerfreiheit der Informationen in diesem Dokument und übernimmt keine Haftung für Fehler, Versäumnisse, Schäden oder Verluste, die aus dem Gebrauch dieses Dokuments entstehen, selbst wenn die Informationen in diesem Dokument genau befolgt werden.

Dieses Dokument ist nicht Teil eines Kaufvertrags zwischen Thermo Fisher Scientific Inc. und einem Kunden. Dieses Dokument regelt oder ändert keine Geschäftsbedingungen; bei widersprüchlichen Informationen zwischen den beiden Dokumenten gelten die Geschäftsbedingungen.

Nur für Forschungszwecke. Dieses Gerät oder Zubehör ist kein Medizinprodukt und ist nicht für die Prävention, Diagnose, Behandlung oder Heilung von Krankheiten vorgesehen.

WARNUNG



Explosions- und Brandgefahren sind zu vermeiden.

Dieses Gerät oder Zubehör ist nicht für den Gebrauch in einer explosionsfähigen Atmosphäre vorgesehen.

Inhalt

| Willkommen | 1 |
|----------------------------|---|
| Typografische Konventionen | 2 |
| Bestellung von Teilen | 3 |
| Kontakt | 3 |

| Erste Schritte mit Ihrem Nicolet Summit OA Spektrometer | 5 |
|---|----|
| Vor dem Eintreffen Ihres Summit Spektrometers | 6 |
| Persönliche Sicherheit mit Ihrem Summit Spektrometer | 9 |
| Auspacken und Einrichten Ihres Spektrometers | 15 |
| Sichern und Wiederherstellen Ihrer Daten | 20 |
| Installieren eines USB-Wireless-Adapters | 27 |
| Anzeigen auf der Lichtleiste des Spektrometers | |
| Anzeigen Ihrer Daten mit der App "OMNIC Anywhere" | |
| | |

| Optionen und Zubehör | 33 |
|---|----|
| Installieren und Warten einer Spüleinheit | 34 |

| Wartung | 45 |
|---|----|
| Wartungsplan für das Summit Spektrometer | 46 |
| Justieren der Spektrometeroptik | 48 |
| Kalibrieren der Laserfrequenz | 49 |
| Reinigen des Spektrometers und Kristalls | 50 |
| Entfernen und Wiedereinsetzen des Kristallaufsatzes | 52 |
| Anbringen der Abdeckung | 54 |
| Auswechseln des Trockenmittels | 59 |
| Austauschen der IR Lichtquelle | 66 |
| | |

Kapitel 1

Willkommen

Mit dem Thermo Scientific Nicolet Summit OA Fourier-Transform-Infrarotspektrometer (FTIR) können chemische Analysen verschiedener Probenmaterialien durchgeführt werden, indem Spektren im mittleren IR-Spektralbereich aufgenommen werden. Das System bietet integrierte Funktionen zur Verifizierung der Messergebnisse, eine leistungsstarke Software-Suite und viele weitere Eigenschaften zur Vereinfachung der Datenaufnahme. Sie können optionale Hardware installieren und verschiedene Service-und Wartungsprozeduren selbst durchführen. Dieses Dokument und die weiterführenden Dokumentationen enthalten alle von Ihnen benötigten Informationen.

Lesen Sie vor der Inbetriebnahme des Spektrometers auch den Standort- und Sicherheitsleitfaden für das Summit Spektrometer.

Hinweis Wir empfehlen, dass Sie die Abdichtung und Trocknung des Spektrometers aufrechterhalten und/oder das Gerät stets spülen. Geräteschäden, die durch die Nichtbeachtung der Maßnahmen zur Abdichtung und Trocknung oder das Unterlassen einer Spülung mit Trockenluft oder Stickstoff entstehen, werden von der Garantie nicht abgedeckt. Wenn Sie Fragen hierzu haben, setzen Sie sich bitte mit uns in Verbindung.

Typografische Konventionen

Sicherheitshinweise und andere wichtige Informationen sind in diesem Handbuch wie folgt gekennzeichnet:

GEFAHR



Gefahren sind zu vermeiden. Weist auf potenziell gefährliche Situationen hin, die bei Nichtvermeidung zu schweren oder tödlichen Verletzungen führen.

WARNUNG



Gefahren sind zu vermeiden. Weist auf potenziell gefährliche Situationen hin, die bei Nichtvermeidung zu schweren oder tödlichen Verletzungen führen können.

VORSICHT



Gefahren sind zu vermeiden. Weist auf eine Gefahrensituation hin, die zu leichten oder mittelschweren Verletzungen führen kann, wenn diese nicht vermieden wird.

HINWEIS

Befolgen Sie die Anweisungen mit dieser Kennzeichnung, um eine Beschädigung des Systems oder einen Datenverlust zu vermeiden.

Hinweis Enthält hilfreiche Zusatzinformationen.

Bestellung von Teilen

Wenden Sie sich für die Bestellung von Teilen an uns.

Wenn Sie das Spektrometer oder ein Zubehör zur Reparatur einsenden möchten, wenden Sie sich bitte zunächst telefonisch oder per E-Mail an uns, damit wir Ihnen die Versanddetails mitteilen oder andere Anweisungen geben können.

Kontakt

Wenden Sie sich für einen technischen Support an www.thermofisher.com.

[Leerseite]

Erste Schritte mit Ihrem Nicolet Summit OA Spektrometer

Bevor Sie die ersten Schritte mit Ihrem neuen Nicolet Summit Lite OA Spektrometer durchführen, sollten Sie vor dem Eintreffen des Geräts die Anforderungen an den Aufstellort sowie die Sicherheit überprüfen, das Gerät auspacken und einrichten sowie jegliche Zubehöre anschließen.

| Vor dem Eintreffen Ihres Summit Spektrometers | |
|--|----|
| Persönliche Sicherheit mit Ihrem Summit Spektrometer | 9 |
| Auspacken und Einrichten Ihres Spektrometers | 15 |
| Sichern und Wiederherstellen Ihrer Daten | 20 |
| Installieren eines USB-Wireless-Adapters | 27 |
| Anzeigen auf der Lichtleiste des Spektrometers | |
| Anzeigen Ihrer Daten mit der App "OMNIC Anywhere" | |

Vor dem Eintreffen Ihres Summit Spektrometers

Bereiten Sie vor dem Eintreffen Ihres Spektrometers einen geeigneten Arbeitsplatz vor. So stellen Sie sicher, dass Sie präzise Messungen durchführen können und das Gerät auch langfristig eine optimale Leistung erreicht.

Dieser Artikel erklärt, welche Schritte Sie durchführen müssen, wenn Ihr Spektrometer eintrifft, und welche Einflüsse (z.B. elektrische Felder oder Umgebungseinflüsse) sich auf die Leistung Ihres Spektrometers auswirken können. Eine ausführliche Erklärung dieser Anforderungen finden Sie im Standort- und Sicherheitsleitfaden für das Spektrometer.

Auspacken des Spektrometers

Wenn Ihr Spektrometer angekommen ist, sollten Sie zwei wichtige Schritte durchführen, bevor Sie den Versandkarton öffnen:

• Untersuchen Sie den Versandkarton auf sichtbare Beschädigungen.

Wenn eine Beschädigung sichtbar ist, wenden Sie sich an uns oder Ihren Händler vor Ort, um die weitere Vorgehensweise zu besprechen.

• Lassen Sie das Spektrometer zunächst im verschlossenen Versandkarton stehen, bis dieses Raumtemperatur angenommen hat.

Das Spektrometer ist im Versandkarton in einer Kunststofftüte verpackt, um es vor Feuchtigkeit zu schützen. Lassen Sie das Spektrometer vor dem Öffnen der Kunststofftüte 24 Stunden stehen, damit es Raumtemperatur annehmen kann. Wenn das Spektrometer ausgepackt wird, bevor es die Raumtemperatur angenommen hat, könnte sich Feuchtigkeit auf den optischen Bauteilen niederschlagen und diese permanent beschädigen.

Schäden, die durch unsachgemäßen Transport oder frühzeitiges Entfernen des Gerätes aus der Kunststofftüte (bevor das Spektrometer Raumtemperatur angenommen hat) verursacht werden, werden nicht durch die Garantie abgedeckt.

Vorbereiten des Arbeitsbereichs

Stellen Sie vor dem Aufstellen des Spektrometers sicher, dass der vorgesehene Arbeitsbereich für das Gewicht und die Abmessungen des Spektrometers geeignet ist. Neben den räumlichen Anforderungen des Spektrometers müssen auch verschiedene Umgebungs- und elektrische Anforderungen berücksichtigt werden.

Abmessungen des Spektrometers

Das Summit Spektrometer benötigt nur eine relativ geringe Stellfläche. Es muss jedoch ausreichend Platz um das Gerät herum vorhanden sein, damit die Wärme aus den Lüftungsschlitzen abgeführt werden kann und die Anschlüsse, der Netzschalter und die Kabel leicht zugänglich sind.

- Gewicht des Summit Spektrometers: 12,6 kg (27,8 lbs)
- Abmessungen (B x H x T): 34 x 24 x 32 cm; (13,3 x 9,6 x 12,7 Zoll)

Umgebungsfaktoren

Das Summit Spektrometer ist ein robustes Gerät, das in verschiedenen Umgebungen eingesetzt werden kann. Die beste Leistung wird jedoch in einer relativ staubfreien Umgebung mit nur geringer Luftfeuchtigkeit erzielt. Das Spektrometer arbeitet zuverlässig in einem Temperaturbereich zwischen 15 °C und 35 °C. Die optimale Leistung wird bei Temperaturen zwischen 20 °C und 22 °C erreicht.

Feuchtigkeit kann innerhalb des Gerätes kondensieren und so die internen Bauteile des Spektrometers beschädigen. Durch das Einhalten einige Vorsichtsmaßnahmen können Sie Ihr Gerät vor Feuchtigkeit schützen:

- Überprüfen Sie den Zustand des Trockenmittels im Gerät in regelmäßigen Abständen, auch wenn dieses über einen längeren Zeitraum nicht verwendet wird.
- Setzen Sie das Gerät keinen raschen Temperaturänderungen aus.
 - Stellen Sie das Gerät nicht in der Nähe von Orten mit heißer oder kalter Luft auf, beispielsweise in die Nähe von Heizungen, Klimaanlagen oder großen Fenstern.

Wenn Ihr Spektrometer in einer relativ feuchten Umgebung aufgestellt werden muss, sollten Sie die Installation eines Spülgasanschlusses in Betracht ziehen.

Elektrische Anforderungen

Die Spannungsversorgung des Spektrometers muss über zweckbestimmte, dauerhaft zur Verfügung stehende Quellen erfolgen, die keine Netzstörungen wie die Folgenden aufweisen:

- Spannungsabfälle
- Spannungsspitzen
- Frequenzschwankungen
- Andere Netzstörungen

Wenn Sie Probleme mit Ihrer Spannungsversorgung vermuten, empfehlen wir eine Überprüfung der Netzqualität. Weitere Informationen erhalten Sie von uns oder von Ihrem Energieversorger vor Ort.

Technische Daten der Spannungsversorgung

Die folgende Tabelle enthält die technischen Daten für die Spannungsversorgung. Sollten Sie Fragen zu den Anforderungen haben, wenden Sie sich an Ihren Service Ansprechpartner vor Ort.

| Anforderungen | Spezifikation |
|--------------------------|--|
| Stromaufnahme | 1,5 A max. |
| Eingangsspannung | AC 100 bis 240 V |
| Netzfrequenz | 50 bis 60 |
| Netzstörungen | Spannungsabfälle und -spitzen sowie andere Störungen dürfen 10 % der Eingangsspannung nicht übersteigen |
| Elektrisches Rauschen | Weniger als 2 V (Gleichtakt) Weniger als 20 V (Gegentakt-Störungen) |

Persönliche Sicherheit mit Ihrem Summit Spektrometer

Obwohl das Nicolet Summit OA Spektrometer als sicheres Spektrometer konstruiert wurde, müssen Sie einige Vorsichtsmaßnahmen beachten, um sich vor potentiellen Gefahren zu schützen, die im täglichen Betrieb und während der Wartung des Gerätes auftreten können.

VORSICHT



Dieses Kapitel ist eine Einführung in die potentiellen Gefahren, denen Sie bei der Nutzung und Wartung des Spektrometers ausgesetzt sein können. Es stellt jedoch keinen umfassenden Leitfaden dar. Bevor Sie das Spektrometer verwenden, sollten Sie die umfassende Beschreibung dieser potentiellen Gefahren im Standort- und Sicherheitsleitfaden für das Spektrometer nachlesen.

Potentielle Gefahren im Normalbetrieb

Während des täglichen Betriebs stellen die folgenden Quellen die größten Gefahren dar:

- Potentiell gefährliche Proben und Lösungsmittel
- Exposition gegenüber Zinkselenid (ZnSe)
- Heiße oder Hitze abstrahlende Teile des Spektrometers
- Der Laser Spektrometers

Sie können Verletzungen oder Beschädigungen des Spektrometers vermeiden, indem Sie die potentiellen Gefahren erkennen und entsprechende Vorsichtsmaßnahmen treffen.

Gefährliche Proben und Lösungsmittel

Treffen Sie entsprechende Vorsichtsmaßnahmen, wenn Sie potentiell gefährliche Proben oder Lösungsmittel verwenden oder analysieren, beispielsweise unter Druck stehende Gase oder korrosive oder brennbare Lösungsmittel.

GEEIGNETE BELÜFTUNG

Es bestehen keine besonderen Anforderungen an die Belüftung Ihres Spektrometers, aber bei der Analyse von potentiell gefährlichen Stoffen kann eine zusätzliche Belüftung erforderlich sein. Stellen Sie sicher, dass bei der Analyse hochgiftiger Proben oder dem Einsatz von Lösungsmitteln, die mit der Infrarot Lichtquelle reagieren oder der Aufnahme von Spektren bei entzündlichen Gasen besondere Belüftungsmaßnahmen getroffen werden.

Bei der Pyrolyse von Lösungsmitteln, die halogenierte Kohlenwasserstoffe enthalten, kann Salzsäure (HCI), Flusssäure (HF) oder Phosgen (COCI₂) entstehen.



FLÜCHTIGE UND ENTFLAMMBARE LÖSUNGSMITTEL

Die Infrarot Lichtquelle im Spektrometer kann entflammbare und flüchtige Proben und Lösungsmittel entzünden. Beachten Sie daher beim Arbeiten mit entflammbaren Proben und Lösungsmitteln die folgenden Sicherheitsmaßnahmen:

- Arbeiten Sie mit eingesetzten Probenraumfenstern.
- Stellen Sie sicher, dass der Arbeitsbereich ordnungsgemäß belüftet wird. Verwenden Sie zu diesem Zweck ein aktives Belüftungssystem ohne Funkenbildung und weitere Zündquellen, welches die Ansammlung entzündlicher Dämpfe in der Luft um das Spektrometer herum unterbindet.
- Bewahren Sie keine entflammbaren Lösungsmittel oder Proben in der Nähe des Spektrometers auf.
- Lassen Sie entflammbare Lösungsmittel oder Proben nicht länger als erforderlich im Probenraum stehen.
- Spülen Sie das Spektrometer mit sauberer Trockenluft oder Stickstoff.

KORROSIVE LÖSUNGSMITTEL

Das Verwenden von Lösungsmitteln im Probenraum, welche HCI- oder HF-Dämpfe entwickeln können, kann das System schwer beschädigen. Wenn Sie halogenierte Lösungsmittel verwenden, spülen Sie das Spektrometer mit sauberer Trockenluft oder Stickstoff.

HINWEIS

Geräteschäden, die durch das Unterlassen einer Spülung oder eine fehlerhafte Spülung des Spektrometers mit Trockenluft oder Stickstoff entstehen, sind von der Garantie ausgenommen.

BIOLOGISCH GEFÄHRLICHE, RADIOAKTIVE UND INFEKTIÖSE SUBSTANZEN

Biologische Proben, z. B. Gewebe, Körperflüssigkeit, infektiöse Substanzen und Blut können Infektionskrankheiten übertragen. Befolgen Sie die Vorgaben Ihrer Organisation zum Arbeiten mit potenziell infektiösen Materialien.

Giftige Materialien

Das Nicolet Summit OA Spektrometer umfasst ein ATR-Kristall aus Zinkselenid (ZnSe), und andere Summit Geräte nutzen u. U. die optionalen ZnSe-Probenraumfenster.

WARNUNG



Das Einatmen und Verschlucken von Giftstoffen ist zu vermeiden.

Zinkselenid (ZnSe) ist giftig. Informationen zur Handhabung und zu Expositionskontrollen entnehmen Sie bitte dem ZnSe-Sicherheitsdatenblatt des Herstellers unter www.specac.com.

Hitzequellen

Teile des Spektrometers können während des täglichen Betriebs sehr heiß werden. Seien Sie insbesondere im Bereich der Infrarot Lichtquelle und den Lüftungsschlitzen vorsichtig.

Die außen liegende Oberfläche der Infrarot Lichtquelle befindet sich an der Unterseite des Spektrometers. Sie kann im Betrieb sehr heiß werden. Die Unterseite des Messgerätes darf während des Betriebs und kurz nach dem Ausschalten nicht berührt werden.

Die Lüftungsschlitze des Spektrometers befinden sich an der linken Seite des Gerätes. Während des normslen Betriebs führt das Spektrometer heiße Luft von den inneren Bauteilen über diese Lüftungsschlitze an die Umgebung ab. Lassen Sie ausreichend Platz um das Spektrometer frei, damit diese heiße Luft abgeführt werden kann.

Laser und optische Sicherheit

Während des normalen Betriebs des Spektrometers sind Sie niemals einer gefährlichen Laserstrahlung ausgesetzt. Wird die Abdeckung des Messgerätes im Rahmen von Wartungsarbeiten entfernt, müssen entsprechende Vorsichtsmaßnahmen getroffen werden, beispielsweise das Tragen einer Schutzbrille. Ihr Servicemitarbeiter wird Sie informieren, wenn dies erforderlich ist.

WARNUNG



Verletzungen sind zu vermeiden.

Blicken Sie niemals direkt in den Laserstrahl oder in reflektierte Laserstrahlung. Nehmen Sie keine Veränderungen am Laser vor und gehen Sie auch beim Austauschen eines defekten Lasers vorsichtig vor, denn Sie könnten dem Laserlicht oder einer hohen Spannung ausgesetzt sein.

Potentielle Gefahren während der Wartung

Im Vergleich zum normalen Betrieb können Sie bei Wartungsarbeiten am Messgerät anderen potentiellen Gefahren ausgesetzt sein. Während der Wartung zählen das Spülen des Messgerätes und das Arbeiten an den internen Bauteilen zu den größten Gefahren.

Spülen des Spektrometers

Insbesondere in feuchten Umgebungen empfehlen wir die Installation einer Spülung für das Spektrometer mit sauberer Trockenluft oder Stickstoff Das Spülen des Spektrometers kann dabei helfen, die internen optischen Bauteile vor Beschädigungen zu schützen, die durch eine feuchte Umgebung oder korrosive Lösungsmittel entstehen. Darüber hinaus hilft es dabei, präzise Ergebnisse sicherzustellen.

Weitere Informationen zum Erwerb und zur Installation einer Spüleinheit für Ihr Spektrometer finden Sie unter "Installieren und Warten einer Spüleinheit".

GEFAHR Vermeiden Sie Feuer und elektrische Schläge.



- Das Spektrometer darf nur mit Stickstoff oder Trockenluft gespült werden.
- Verwenden Sie zum Spülen dieses Geräts niemals entzündliche, brennbare oder toxische Gase. Das Spülgas muss frei von Öl und anderen reaktiven Materialien sein. Durch die von der Lichtquelle ausgehende oder durch Laserabsorption entstehende Wärme können brennbare Gase oder reaktive Materialien im Spülgas entzündet werden.

Arbeiten an internen Bauteilen

In der Regel gibt es keinen Grund, die Abdeckung des Spektrometers zu entfernen oder zu versuchen, an die internen Bauteile zu gelangen. Wenn Sie die Abdeckung jedoch für Wartungsarbeiten entfernen müssen, besteht die Gefahr eines (elektrischen) Schocks, Verbrennungen oder der Aussetzung gegenüber Laserlicht. verursachen.

VORSICHT



VORSICHT



Brandgefahren sind zu vermeiden.

Verletzungen sind zu vermeiden.

Gefahren durch Stromschläge sind zu vermeiden.

Interne Bauteile, insbesondere die Infrarot Lichtquelle, können während des normalen Betriebs sehr heiß werden. Schalten Sie das Messgerät aus und warten Sie mindestens 10 Minuten, bevor Sie Bauteile auswechseln.

Nachdem das Gerät von allen Spannungsquellen getrennt wurde, können die Kondensatoren bis zu 30 Sekunden lang geladen bleiben und einen Stromschlag

WARNUNG



- Blicken Sie niemals direkt in den Laserstrahl oder in reflektierte Laserstrahlung. Veränderungen am Laser sind nicht zulässig. Sie könnten dem Laserlicht oder hoher Spannung ausgesetzt werden.
- Wenn Sie den Laser justieren oder Verfahren durchführen, die nicht in der Bedienungsanleitung oder den Handbüchern beschrieben sind, können Sie einer gefährlichen Strahlung ausgesetzt werden.

AUSWECHSELN DES TROCKENMITTELS

Beim Öffnen des Trockenmittelfachs müssen Sie verhindern, dass entflammbare Flüssigkeiten oder Gase in das Fach eintreten. Anweisungen zum Auswechseln des Trockenmittels finden Sie unter "Auswechseln des Trockenmittels".

GEFAHR

Die Gefahr einer Explosion ist zu vermeiden.



Schalten Sie das Messgerät stets aus, ziehen Sie den Netzstecker und entfernen Sie die Proben aus dem Gerät, bevor Sie das Trockenmittelfach öffnen. Das Eindringen von entflammbaren Flüssigkeiten und Gasen in das Trockenmittelfach stellt eine Explosionsgefahr dar. Sollte dieser Fall eintreten, setzen Sie sich umgehend mit uns in Verbindung. Das Gerät darf nicht an die Spannungsversorgung angeschlossen werden, bis das Problem behoben wurde.

AUSTAUSCHEN DER IR LICHTQUELLE

Die Infrarot Lichtquelle wird während des Normalbetriebs sehr heiß. Wenn Sie die IR Quelle austauschen müssen, schützen Sie sich vor Verbrennungen und beachten Sie die Explosionsgefahr.

VORSICHT



Brandgefahren sind zu vermeiden.

Die IR Quelle wird während des Normalbetriebs des Spektrometers sehr heiß. Lassen Sie die IR-Quelle nach dem Ausschalten mindestens 10 Minuten lang abkühlen, bevor Sie Arbeiten daran vornehmen.

GEFAHR



Die Gefahr einer Explosion ist zu vermeiden.

Bevor Sie die IR Quelle aus dem Spektrometer entfernen, müssen Sie das Spektrometer ausschalten, den Netzstecker ziehen und die Proben aus dem Gerät entfernen. Das Eintreten von entflammbaren Flüssigkeiten und Gasen in das Fach der IR Quelle stellt eine Explosionsgefahr dar. Sollte dieser Fall eintreten, setzen Sie sich umgehend mit uns in Verbindung. Das Gerät darf nicht an die Spannungsversorgung angeschlossen werden, bis das Problem behoben wurde.

Zusammenfassung

Das Nicolet Summit OA Spektrometer ist ein sicheres und robustes Messgerät, dennoch bestehen während des normalen Betriebs und der Wartung bestimmte Gefahren. Während des normalen Betriebs muss mit potentiell gefährlichen Proben und Lösungsmitteln vorsichtig vorgegangen werden, und die heißen oder Hitze abstrahlenden Teile des Spektrometers sind zu meiden. Während der Wartung sind entsprechende Vorsichtsmaßnahmen zu treffen, um Verletzungen oder Beschädigungen beim Spülen des Messgerätes, beim Handhaben der internen Bauteile und beim Auswechseln des Trockenmittels zu vermeiden.

Auspacken und Einrichten Ihres Spektrometers

Das Auspacken und Einrichten Ihres Thermo Scientific Nicolet Summit OA Spektrometers umfasst das Aufstellen des Spektrometers am vorgesehenen Arbeitsbereich, das Anschließen und Einschalten des Spektrometers und das Durchführen eines kurzen Leistungstests.

Bevor Sie beginnen

HINWEIS

Warten Sie 24 Stunden, bevor Sie das Spektrometer aus der Plastiktüte nehmen, um zu verhindern, dass sich Feuchtigkeit auf der internen Optik niederschlägt und diese beschädigt. Schäden, die aufgrund eines frühzeitigen Öffnen der Plastiktüte verursacht werden (bevor das Spektrometer die Raumtemperatur angenommen hat), sind nicht durch die Garantie abgedeckt.

Zur Erinnerung: Lesen Sie den Standort- und Sicherheitsleitfaden für das Summit Spektrometer, in dem Sie weitere Informationen zu einem sicheren und geeigneten Aufstellort für Ihr Messgerät finden.

Auspacken und Einrichten

Die Schritte zum Auspacken und Einrichten Ihres Spektrometers hängen von den Optionen ab, die Sie mit dem Gerät erworben haben.

1. Auspacken des Versandkartons

Entnehmen Sie das Spektrometer aus dem Versandkarton und entfernen Sie die zusätzlichen Verpackungsmaterialien.

Das Spektrometer ist mit einer abnehmbaren Staubabdeckung ausgestattet, die über dem ATR-Kristall sitzt. Verwenden Sie die Staubabdeckung bei der Lagerung und beim Transport des Spektrometers und entfernen Sie sie für den regulären Gebrauch.

- 2. Schließen Sie das Spektrometer an eine Spannungsversorgung und an eine ordnungsgemäß geerdete Wechselspannungsquelle an.
- 3. Schließen Sie Monitor, Tastatur und Maus an.

- a. Schließen Sie den Monitor an den Mini DisplayPort-Anschluss des Spektrometers an.
- b. Schließen Sie die Tastatur und die Maus an die USB-Anschlüsse des Spektrometers an.

Hinweis Einige USB- und Mini DisplayPort-Geräte sind bekannt dafür, dass sie Signale von drahtlosen USB-Geräten stören, beispielsweise von drahtlosen Tastaturen und Mäusen. Wenn Sie Probleme mit Ihrer drahtlosen Tastatur oder Ihrer drahtlosen Maus bemerken, beispielsweise Verzögerungen oder nicht übertragene Eingabebefehle, verwenden Sie das mitgelieferte USB Verlängerungskabel, um den Empfänger für die drahtlosen USB Geräte weiter entfernt von den USB Anschlüssen zu positionieren.

4. Drücken Sie den Netzschalter am Spektrometer, um das Gerät einzuschalten. Die Lichtleiste des Spektrometers zeigt ein grünes Lauflicht, während das Gerät aufheizt. Ist das Spektrometer betriebsbereit, wird die Lichtleiste als ein durchgehend grüner Balken angezeigt.

Beim Ein- oder Ausschalten des Spektrometers hören Sie u. U. das Geräusch, das beim Bewegen des internen Polystyrol-Vergleichsstandards entsteht. Dies ist normal.



Wenn Sie das Spektrometer zum ersten Mal einschalten, werden Sie aufgefordert ein neues Kennwort einzugeben. Wir empfehlen, dass Sie zu diesem Zeitpunkt kein Kennwort eingeben. Fahren Sie zunächst ohne Eingabe eines Kennworts fort und kümmern Sie sich dann später, nachdem Sie die Einrichtung abgeschlossen haben, um die Verwaltung der Konten und Kennwörter unter Berücksichtigung der Richtlinien Ihrer Organisation.

Hinweis Wenn der Bildschirm nach dem Einschalten des Systems schwarz bleibt, überprüfen Sie die Stromversorgung des Monitors, den DisplayPort-Anschluss und die Monitoreinstellungen. Weitere Informationen zum Ändern der Einstellungen finden Sie in der Dokumentation Ihres Monitors. 5. Öffnen Sie die OMNIC Paradigm-Software.

Es kann mehrere Minuten dauern, bis die Software gestartet ist und eine Verbindung zum Spektrometer hergestellt hat. Der Verbindungsstatus wird oben rechts auf dem Bildschirm angezeigt. Wenn der Status anzeigt, dass eine Verbindung hergestellt wurde, fahren Sie mit dem nächsten Schritt fort.



Hinweis Wenn Sie vor der Herstellung einer Verbindung zum Spektrometer Menüs in der Software öffnen, werden die Justierung und Kalibrierung deaktiviert. Um diese Funktionen zu aktivieren, klicken Sie in der Symbolleiste auf Dashboard, wenn eine Verbindung zum Spektrometer hergestellt ist.

- 6. Überprüfen der Spektrometerleistung
 - a. Justieren Sie die Spektrometeroptik.
 - i. Wählen Sie auf der Desktop-Bedienungsoberfläche Daten erfassen > Diagnostik
 > Spektrometer justieren.
 - ii. Wählen Sie **Start**, um mit der Justierung zu beginnen. Bei erfolgreichem Abschluss der Justierung wird auf dem Bildschirm "Justierung abgeschlossen" angezeigt.
 - b. Kalibrieren Sie den Laser.
 - i. Wählen Sie Daten erfassen > Diagnostik > Laserkalibrierung.
 - ii. Wählen Sie Start, um mit der Kalibrierung zu beginnen.
 - iii. Die Laserkalibrierung nimmt einige Minuten in Anspruch. Nach Abschluss der Kalibrierung zeigt das Display die Meldung "Messung abgeschlossen" an und meldet ggf. Änderungen in der Laserfrequenz.

Das Spektrometer braucht bis zu 12 Stunden, um ein thermisches Gleichgewicht zu erreichen. Um die bestmöglichen Ergebnisse zu erzielen, führen Sie die Ausrichtung und Kalibrierung nach den ersten 12 Betriebsstunden erneut aus.

c. Führen Sie den Arbeitsablauf Nicolet Summit OA - PV-Test aus.

Im Rahmen des Arbeitsablaufs der Leistungsüberprüfung wird ein Referenzmaterial verwendet und eine Reihe von Standardtests durchgeführt, um die Betriebsbereitschaft des Spektrometers sowie die Genauigkeit der Messdaten zu überprüfen.

- i. Führen Sie vom Dashboard der Desktop-Bedienungsoberfläche einen Bildlauf zum Bereich "Arbeitsabläufe" durch.
- ii. Klicken Sie im Bereich "Arbeitsablauf" mit der rechten Maustaste auf den Arbeitsablauf Nicolet Summit OA PV und wählen Sie "Ausführen".



- iii. Folgen Sie den Anweisungen auf dem Bildschirm, um den Arbeitsablauf abzuschließen.
- iv. Nach der Anzeige der Tabelle "PV-Ergebnisse" wählen Sie "Fertig stellen", um den Arbeitsablauf abzuschließen und zum Dashboard zurückzukehren, oder wählen Sie "Squalan-Spektrum messen", um mit dem optionalen Squalan-Teil des Arbeitsablaufs fortzufahren. Die Messung eines Squalan-Spektrums ist eine weitere Vergleichsprüfung, um die Leistung des Systems zu überprüfen.



Nächste Schritte

Herzlichen Glückwunsch! Ihr Spektrometer ist jetzt einsatzbereit. Bevor Sie mit der Aufnahme von Spektren beginnen, sollten Sie eine Sicherheitskopie Ihres Systems anlegen und ein

Erste Schritte mit Ihrem Nicolet Summit OA Spektrometer

Wiederherstellungslaufwerk erstellen, um Ihr System vor möglichen Datenverlusten in der Zukunft zu schützen. Windows 10 umfasst verschiedene integrierte Tools zur Sicherung und Wiederherstellung Ihrer Daten.

Weitere Informationen zum Erstellen eines Wiederherstellungslaufwerks und Anlegen eines Systemabbilds sowie zum Wiederherstellen Ihres Systems nach einem Datenverlust finden Sie unter "Sichern und Wiederherstellen Ihrer Daten".

Weitere Informationen zur Identifizierung und Analyse von Proben mit dem Spektrometer finden Sie in den OMNIC Paradigm-Bedienungsanleitungen und -Tutorials unter thermofisher.com/ftir-help.

Zum Wechseln zur Bedienoberfläche wählen Sie Ansicht / Anzeige > Bediener.

Detaillierte Informationen zum Erstellen und Umsetzen individueller Lösungen mit dem Package Editor und der Bedienoberfläche finden Sie unter "Individuelle Lösungen" in der Bedienungsanleitung der OMNIC Paradigm Software.

Sichern und Wiederherstellen Ihrer Daten

Bevor Sie mit der Datenaufnahme beginnen, sollten Sie einige Vorsichtsmaßnahmen treffen, um Ihr System vor möglichen Datenverlusten in der Zukunft zu schützen. Windows 10 umfasst verschiedene integrierte Tools zur Sicherung und Wiederherstellung Ihrer Daten.

WICHTIG! Um sicherzustellen, dass Sie im Falle einer Beschädigung der Festplatte oder eines Ausfalls des Betriebssystems alle Daten wiederherstellen können, sollten Sie Wiederherstellungsmedien erstellen, ein Systemabbild speichern und ein Verfahren zur Sicherung Ihrer Daten festlegen.

Erstellen eines Wiederherstellungslaufwerks

Erstellen Sie ein Wiederherstellungslaufwerk, damit Sie bei Problemen mit Ihrem Betriebssystem auf die erweiterten Startoptionen von Windows 10 zugreifen können - z. B. Wiederherstellung eines Systemabbilds, Systemwiederherstellung und Starteinstellungen - selbst wenn Ihre Windows-Installation beschädigt wurde.

Erforderliche Materialien

• Ein FAT32 formatiertes Flashlaufwerk mit mindestens 8 GB Speicherplatz

HINWEIS

Alle auf diesem Laufwerk vorhandenen Daten werden während dieses Prozesses gelöscht. Stellen Sie, bevor Sie fortfahren, sicher, dass Sie eine Sicherheitskopie aller Daten auf dem Flashlaufwerk erstellt haben, die Sie behalten möchten.

So erstellen Sie ein Wiederherstellungslaufwerk:

- 1. Schließen Sie das Flashlaufwerk an einen der USB Anschlüsse an der Seite des Spektrometers an.
- Geben Sie in das Suchfeld in der Taskleiste von Windows 10 den Begriff "Wiederherstellungslaufwerk erstellen" ein und wählen Sie Wiederherstellungslaufwerk erstellen.

Erste Schritte mit Ihrem Nicolet Summit OA Spektrometer



- 3. Wenn Sie gefragt werden, ob Sie Änderungen durch diese App zulassen möchten, wählen Sie Ja.
- 4. Deaktivieren Sie das Kontrollkästchen "Sichert die Systemdateien auf dem Wiederherstellungslaufwerk" und klicken Sie auf **Weiter**.
- 5. Wählen Sie das angeschlossene Flashlaufwerk in der Liste der verfügbaren Laufwerke aus und klicken Sie auf **Weiter**.



- 6. Klicken Sie auf **Erstellen**. Wenn der Vorgang abgeschlossen ist, klicken Sie auf **Fertig stellen**, um das Fenster zu schließen.
- 7. Entfernen Sie das Flashlaufwerk und bewahren Sie es an einem sicheren Ort auf. Sie sollten das Laufwerk mit einem aussagekräftigen Namen wie "Summit-Wiederherstellungslaufwerk" beschriften.

Weitere Informationen zum Erstellen von Wiederherstellungsmedien finden Sie in der Online-Hilfe von Microsoft.

Anlegen eines Systemabbilds

Ein Systemabbild ist eine Momentaufnahme Ihres gesamten Systems, einschließlich Ihrer Daten und Dateien, installierter Anwendungen und anderer Einstellungen. Bei einem vollständigen Ausfall der Festplatte oder des Betriebssystems können Sie Ihr System mit dem Systemabbild wieder in den Zustand versetzen, in dem es beim Erstellen des Abbildes war. Es liegt an Ihnen, wie oft Sie das Systemabbild erstellen bzw. aktualisieren. Wir empfehlen, zumindest ein Abbild einer sauberen Installation nach der Ersteinrichtung des Spektrometers anzulegen.

Je nachdem, wie viele Daten Sie in das Systemabbild kopieren, kann dieser Vorgang eine Weile dauern. Wenn Sie eine große Datenmenge kopieren, kann dieser Vorgang mehr als eine Stunde in Anspruch nehmen.

Erforderliche Materialien

- NTFS-formatierte SSD
- So erstellen Sie ein Systemabbild:
- 1. Schließen Sie die SSD an einen der USB Anschlüsse des Spektrometers an.
- Öffnen Sie in Windows 10 die Systemsteuerung und wählen Sie Sichern und Wiederherstellen (Windows 7).
- 3. Wählen Sie im linken Bereich die Option Systemabbild erstellen aus.

| 🛞 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. | | |
| | Select another backup to restore files from | | |
| | | | |
| | | | |
| | | | |

4. Wählen Sie Auf Festplatte und wählen Sie Ihre SSD in der Liste aus. Klicken Sie auf Weiter.

Erste Schritte mit Ihrem Nicolet Summit OA Spektrometer

| 虪 Create a system image | |
|---|---|
| Where do you want to save the backu | p? |
| A system image is a copy of the drives required f drives. A system image can be used to restore yo stops working; however, you can't choose indivic | r Windows to run. It can also include additional r computer if your hard drive or computer ever ual items to restore. |
| On a hard disk | |
| - Backup (E:) 465.60 GB free | ~ |
| | |
| ○ On one or more DVDs | |
| ○ On one or more DVDs | |
| On one or more DVDs DVD Drive (D:) On a network location | v |
| 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 (D:) On a network location | ∽ Select |
| On one or more DVDs DVD Drive (D:) On a network location | ✓ Select |
| O On one or more DVDs DVD Drive (D:) On a network location | ∽ Select |

- 5. Bestätigen Sie die Datensicherungseinstellungen und klicken Sie auf Sicherung starten.
- 6. Klicken Sie bei der Frage "Möchten Sie einen Systemreparaturdatenträger erstellen?" auf Nein.
- Klicken Sie auf Schließen und entfernen Sie die SSD vom Summit Spektrometer. Beschriften Sie die SSD beispielsweise mit "Summit-Systemabbild" und bewahren Sie diese an einem sicheren Ort auf.

Wiederherstellen Ihres Systems

Bei einem Ausfall des Betriebssystems verwenden Sie Ihr Wiederstellungslaufwerk und Ihr Systemabbild, um Ihr System wiederherzustellen.

Erforderliche Materialien

- USB Wiederherstellungslaufwerk
- SSD mit Ihrem Systemabbild

So stellen Sie Ihr System mithilfe Ihres Wiederstellungslaufwerks und Ihres Systemabbilds wieder her:

- 1. Schalten Sie das Summit Spektrometer aus.
- 2. Schließen Sie eine externe Tastatur an einen der USB Anschlüsse des Summit Spektrometers an.
- 3. Schließen Sie das USB Wiederherstellungslaufwerk an einen der USB Anschlüsse an.
- 4. Schalten Sie das System ein.

5. Wenn der Thermo Scientific-Willkommenbildschirm angezeigt wird, drücken Sie mehrmals die **Esc**-Taste auf der Tastatur, um das System-BIOS anzuzeigen.

Wenn das Popupfenster "Ohne Speichern beenden ?" angezeigt wird, wählen Sie mit den Pfeiltasten die Option "Nein".

- 6. Navigieren Sie mithilfe der Pfeiltaste zur Registerkarte Speichern und Beenden.
- 7. Navigieren Sie zum Abschnitt "Boot Reihenfolge ändern" und wählen Sie Ihr Wiederherstellungslaufwerk. Drücken Sie die **Eingabetaste**, um vom Wiederherstellungsabbild zu booten.



- 8. Wenn der Bildschirm "Auswahl des Tastaturlayouts" angezeigt wird, entfernen Sie das Flash Wiederherstellungslaufwerk vom USB Anschluss und schließen die SSD mit dem Systemabbild an.
- 9. Wählen Sie das von Ihnen bevorzugte Tastaturlayout.
- 10. Wählen Sie im Bildschirm "Eine Option auswählen" die Option Fehlerbehebung.



- 11. Wählen Sie Erweiterte Optionen.
- 12. Wählen Sie unter "Erweiterte Optionen" die Option Wiederherstellung eines Systemabbilds.

Erste Schritte mit Ihrem Nicolet Summit OA Spektrometer



- 13. Wählen Sie "Windows 10" als Zielbetriebssystem.
- 14. Wählen Sie im Fenster "Wählen Sie ein Systemabbild aus" das letzte verfügbare oder ein anderes Systemabbild aus. Klicken Sie auf **Weiter**.



 Klicken Sie auf Weiter, um die erweiterten Wiederherstellungsoptionen auf den Standardeinstellungen zu belassen, und klicken Sie auf Fertig stellen, um das Systemabbild wiederherstellen. Klicken Sie zur Bestätigung auf Ja.

Die Wiederherstellung dauert etwa 15 Minuten, und das System wird nach Abschluss des Vorgangs automatisch neu gebootet.

Wenn das System wiederhergestellt wurde, entfernen Sie die SS vom Spektrometer und bewahren das Wiederherstellungslaufwerk und die SSD mit dem Systemabbild wieder an einem sicheren Ort auf.

Weitere Informationen zum Wiederherstellung des Systems aus einem Systemabbild finden Sie in der Online-Hilfe von Microsoft.

Sicherung Ihrer Systemdateien

Zusätzlich zum Erstellen eines Wiederherstellungsmediums und eines Systemabbilds sollten Sie auch Ihre Daten in regelmäßigen Abständen sichern. Windows 10 bietet verschiedene Optionen zum Sichern Ihrer Dateien, beispielsweise das Tool "Dateiversionsverlauf". Das IT-Personal Ihrer Organisation hält weitere Informationen bereit, welche Sicherungsmethode für die Anforderungen Ihrer Organisation am besten geeignet ist.

Installieren eines USB-Wireless-Adapters

Sie können das Nicolet Summit OA Spektrometer mit einem optionalen USB-Wireless-Adapter an ein Wireless-Netzwerk anschließen. Wenn Sie den Adapter mit Ihrem Spektrometer erworben haben, ist er bereits vorinstalliert.

Zum Installieren des USB-Wireless-Adapters schließen Sie den Adapter an einen USB-Anschluss an. Für eine längerfristige Installation schließen Sie den Adapter an einen Anschluss unter der Geräteabdeckung an.

So installieren Sie einen USB-Wireless-Adapter:



1. Schließen Sie den Adapter an einen verfügbaren USB-Anschluss an.

2. Der Adapter wird vom System automatisch erkannt. Informationen zum Verbinden des Adapters mit einem Netzwerk finden Sie im Support von Microsoft.

So schließen Sie den Adapter an einen USB-Anschluss unter der Geräteabdeckung an:

Für die Installation müssen der Kristallaufsatz und die Geräteabdeckung entfernt werden.

- 1. Anweisungen zum Entfernen des Kristallaufsatzes und der Geräteabdeckung finden Sie unter <u>"Anbringen der Abdeckung"</u>.
- 2. Schließen Sie den Adapter bei abgenommener Abdeckung an einen der verfügbaren USB-Anschlüsse an.

Erste Schritte mit Ihrem Nicolet Summit OA Spektrometer



3. Setzen Sie die Geräteabdeckung und den Kristallaufsatz wieder ein.

Das Spektrometer ist nun bereit für den Anschluss an ein Wireless-Netzwerk.

Anzeigen auf der Lichtleiste des Spektrometers

Das Thermo Scientific Nicolet Summit OA Spektrometer umfasst eine leicht ablesbare Lichtleiste (die "LightBar"), auf der Sie mit einem Blick erkennen können, ob ein Produkt die Qualitätsprüfungen bestanden hat. Darüber hinaus zeigt die Lichtleiste den Status des Spektrometers an.

In den folgenden Tabellen sind die verschiedenen Lichtsignale der Lichtleiste und deren Bedeutung aufgeführt.

AUFHEIZEN

| Signal | Systemstatus | Beschreibung |
|------------------|--|---|
| Grünes Lauflicht | Das System ist eingeschaltet und heizt auf | Das grüne Licht bewegt sich kontinuierlich von links nach rechts. Das Aufheizen dauert etwa 2 Minuten. |

HERKÖMMLICHE MESSUNG

| Signal | Systemstatus | Beschreibung |
|----------------|-----------------------------|---|
| Dauerhaft grün | Betriebsbereit | Die gesamte Lichtleiste leuchtet grün |
| Grüner Komet | Daten werden aufgenommen | Ein kleiner Bereich der grünen Lichtleiste bewegt sich hin und her |

SMARTE MESSUNG

| Signal | Systemstatus | Beschreibung |
|--------------|--|---|
| Blauer Puls | Smarte Messung des Backgrounds | Die gesamte Lichtleiste pulsiert blau |
| Blauer Komet | Probenspektrum wird gemessen (vom Anwender initiiert) | Ein kleiner Bereich der blauen Lichtleiste bewegt sich hin und her |

ANALYSEERGEBNISSE

| Signal | Systemstatus | Beschreibung |
|------------------|--|---|
| Grüne %-Anzeige | Gütewert oder QCheck- Ergebnis oberhalb des definierten Schwellwerts | Die Anzahl an grünen Segmente entspricht dem Gütewert. Beispielsweise entsprechen etwa 90 % grüne Segmente einem Gütewert von 90. |
| Orange %-Anzeige | Gütewert oder QCheck- Ergebnis unterhalb des definierten Schwellwerts | Die Anzahl an orangefarbenen Segmente entspricht dem Gütewert. Beispielsweise entsprechen etwa 30 % grüne Segmente einem Gütewert von 30. |

DIAGNOSE

| Signal | Systemstatus | Beschreibung |
|---------------|--------------|---|
| Rotes Blinken | Systemfehler | Die gesamte Lichtleiste blinkt rot. Weitere Informationen zur Beschreibung des Fehlers finden Sie unter Systemstatus in der OMNIC Paradigm Software. |

Anzeigen Ihrer Daten mit der App "OMNIC Anywhere"

OMNIC Anywhere ist eine Cloud-basierte Anwendung, mit der Sie Ihre Daten auf einem angeschlossenen PC, Apple Computer oder Android- bzw. iOS-Gerät anzeigen können.



Abbildung 2-1: OMNIC Anywhere in einem Webbrowser

Mit einem Nicolet Summit OA Spektrometer und einem kostenlosen Connect-Account können Sie Proben in einem Klassenraum oder Labor messen und die aufgenommenen Daten auf Ihren Connect-Account hochladen, um sie dann auf einem anderen Gerät auf dem Campus oder in einer anderen Abteilung anzuzeigen, zu analysieren oder zu teilen.

Weitere Informationen zum Erstellen eines Connect-Accounts und zum Anzeigen Ihrer Daten mit OMNIC Anywhere finden Sie unter <u>Cloud-enabled FTIR Spectroscopy</u>.

Kapitel 3

Optionen und Zubehör

Das Nicolet Summit OA kann mit der optionalen Spüleinheit betrieben werden.

| Installieren und Warten einer | Spüleinheit |
|-------------------------------|-------------|
|-------------------------------|-------------|
Installieren und Warten einer Spüleinheit

Das Spülen des Spektrometers mit Trockenluft oder Stickstoff schützt die internen Bauteile vor Feuchtigkeit und anderen Fremdstoffen aus der Umgebung. Das Installieren einer Spüleinheit erfordert das Montieren von Ventilen und Reglern, das Austauschen der Trockenmittelkartusche, das Anschließen des Spülgases und das Einstellen des Drucks und der Flussrate.

WARNUNG

Explosionsgefahren sind zu vermeiden.



Verwenden Sie zum Spülen dieses Geräts niemals entzündliche, brennbare oder toxische Gase. Das Spülgas muss frei von Öl und anderen reaktiven Materialien sein. Durch die von der Lichtquelle ausgehende oder durch Laserabsorption entstehende Wärme können brennbare Gase oder reaktive Materialien im Spülgas entzündet werden. Das Gerät darf nur mit Stickstoff oder getrockneter Luft gespült werden.

HINWEIS

Wir empfehlen, dass Sie die Vorrichtungen zur Abdichtung und Trocknung aufrechterhalten und/oder das Gerät stets spülen. Die Garantie deckt keine Schäden am Gerät ab, die durch das Unterlassen einer ordnungsgemäßen Trocknung oder Spülung des Spektrometers entstehen. Wenn Sie Fragen hierzu haben, setzen Sie sich bitte mit uns in Verbindung.

ERFORDERLICHES SETUP:

Bevor Sie die Spüleinheit installieren, müssen Sie eine Quelle für Trockenluft oder Stickstoff vorsehen, die den Spezifikationen von Thermo Scientific für das Nicolet[™] Summit Spektrometer entsprechen. Weitere Informationen können Sie dem Standort- und Sicherheitsleitfaden für das Summit Spektrometer entnehmen.

Hinweis Für ein optimales Spülergebnis trocknen Sie das Spülgas auf einen Taupunkt von -70 °C (-94 °F) oder kälter.

ERFORDERLICHE MATERIALIEN:

Zusätzlich zur Spüleinheit benötigen Sie die folgenden Werkzeuge und Materialien:

- 3/4-Zoll-Maulschlüssel
- 11/16-Zoll-Maulschlüssel
- Einen #2 Phillips Kreuzschlitzschraubendreher

- 3/32-Zoll-Innensechskantschlüssel
- 7/64-Zoll-Innensechskantschlüssel
- Gewindedichtband ("Klempnerband" oder "Teflonband")

So installieren Sie eine Spüleinheit:

- 1. Befestigen der Rohrleitungsbaugruppe an der Spülgasquelle
 - a. Bringen Sie den Regler und entweder ein 1/4-Zoll-Fitting mit Außengewinde oder ein 3/8-Zoll-Fitting mit Innengewinde an der Spülgasquelle an. (Wählen Sie den Regler und die Fittings passend zur Spülgasquelle aus.)



b. Wenn Sie ein 1/4-Zoll-Fitting mit Außengewinde verwenden, gehen Sie zum nächsten Schritt.

Wenn Sie ein 3/8-Zoll-Fitting mit Innengewinde an der Spülgasquelle verwenden, montieren Sie das Reduzierstück (3/8 auf 1/4 Zoll), das mit der Spüleinheit geliefert wurde. Umwickeln Sie das Gewinde des Reduzierstücks vor der Montage mit Teflon-Dichtband und ziehen Sie die Verbindung mit einem 11/16-Zoll-Maulschlüssel fest.



c. Umwickeln Sie das Gewinde des Reduzierstücks oder des 1/4-Zoll-Fittings mit Außengewinde mit Teflon-Dichtband und schließen Sie dann die Druckkupplung an. Ziehen Sie die Verbindung mit einem 3/4-Zoll-Maulschlüssel fest.



- d. Lassen Sie den Einlass der Rohrleitungsbaugruppe fest in die Schnellkupplung einrasten.
- 2. Schalten Sie das Spektrometer aus.
 - a. Drücken Sie zum Ausschalten des Spektrometers die Netztaste.
 - b. Ziehen Sie das Netzkabel ab.
- 3. "Entfernen des Kristallaufsatzes"

- 4. Nehmen Sie die Geräteabdeckung ab.
 - a. Neigen Sie das Spektrometer vorsichtig auf seine Rückseite, sodass es auf den Rippen des Kühlkörpers zum Liegen kommt.
 - b. Lösen Sie mit einem 7/64-Innensechskantschlüssel die vier Schrauben, mit denen die Abdeckung an der Grundplatte des Geräts befestigt ist.



c. Stellen Sie das Spektrometer vorsichtig wieder auf die Füße. Die vier Schrauben fallen heraus, wenn Sie das Spektrometer wieder auf die Füße stellen. Achten Sie darauf, dass sie nicht verloren gehen.

d. Heben Sie die Abdeckung in einer geraden Bewegung nach oben ab.



- 5. Entfernen der Trockenmittelkartusche
 - a. Lösen Sie die beiden Schrauben mit dem #2 Philipps Kreuzschlitzschraubenzieher von der Trockenmittelkartusche, und ziehen Sie diese gerade nach oben aus dem Spektrometer heraus.



b. Entfernen und entsorgen Sie den großen O-Ring.



Hinweis Achten Sie darauf, dass der O-Ring nicht in die Trockenmittelkammer fällt.

- 6. Öffnen Sie die Spüleinheit und das versiegelte Trockenmittelpaket.
- Entfernen Sie den O-Ring aus dem Paket und montieren Sie ihn an der Grundplatte des Messgerätes (siehe vorherige Abbildung). Drücken (nicht rollen) Sie den neuen O-Ring in die richtige Position, um sicherzustellen, dass er korrekt in der Nut sitzt.
- 8. Entnehmen Sie die Trockenmittelkartusche mit den Spülanschlüssen aus dem Paket, setzen Sie die Kartusche in das Trockenmittelfach ein und stellen Sie sicher, dass sie korrekt über dem O-Ring sitzt. Dann fixieren Sie die Kartusche mit den beiden Schrauben.

Achten Sie auf eine korrekte Ausrichtung der Kartusche. Das Etikett auf der Kartusche sollte nach dem Einbau der Kartusche von der Vorderseite des Spektrometers aus gelesen werden können.



9. Verlegen Sie die Spülleitung durch die Spülleitungsöffnung auf der Rückseite des Spektrometers.





- b. Senken Sie die Abdeckung vollständig ab. Wenn die Abdeckung vollständig abgesenkt ist, sitzt die Spülleitung in dem Schlitz zwischen Abdeckung und Kühlkörper.
- c. Verlegen Sie die Spülleitung so, dass sie zur Seite verläuft und neigen Sie das Spektrometer

dann auf die Rückseite, sodass es auf den Rippen des Kühlkörpers zum Liegen kommt.

- d. Setzen Sie die vier Schrauben ein, die Sie zuvor entfernt haben. Ziehen Sie alle Schrauben fest.
- e. Stellen Sie das Spektrometer wieder auf die Füße.
- f. Schieben Sie die überschüssige Länge der Leitung in das Spektrometer oder ziehen Sie die Leitung bei Bedarf weiter heraus.



- 10. "Wiedereinsetzen des Kristallaufsatzes".
- 11. Schließen Sie die Durchflusskupplung von der Rohrleitungsbaugruppe (mit der Kennzeichnung "zum Spektrometer") an die Spülleitung an.
- 12. Einstellung der Spülgasregelung
 - a. Öffnen Sie das Hauptreglerventil und stellen Sie die Spülgasregelung wie im Folgenden gezeigt ein (siehe Einstellen der Spülgasregelung).

| Hardware | Druck (psig) | Durchflussrate (scfh) |
|---------------------|--------------|-----------------------|
| Summit Spektrometer | 5 | 1 |

- b. Schließen Sie das Netzkabel an die Netzsteckdose an und schalten Sie das Gerät am Netzschalter ein.
- c. Warten Sie 30-60 Minuten, damit das Spektrometer vollständig gespült ist.

Einstellen der Spülgasregelung

Eine korrekte Einstellung der Spülgasregelung schützt Ihr Spektrometer vor Feuchtigkeit, ohne das Gerät Vibrationen auszusetzen.

Ein optimales Spülergebnis wird mit Spülgas erzielt, das auf einen Taupunkt von 70 °C oder darunter abgekühlt wird.

So stellen Sie die Spülgasregelung ein:

- 1. Öffnen Sie das Hauptventil, um den Fluss des Spülgases durch den Regler zu starten.
- 2. Stellen Sie den Druckregler für das Spektrometer so ein, dass das Manometer einen Druck von 34 kPa (5 psig) anzeigt.
- 3. Stellen Sie den Durchflussmesser für das Spektrometer auf einen Wert von 0,47 l/min (1 scfh) ein.

HINWEIS

Flussraten über den empfohlenen Werten können Vibrationen verursachen, die die Qualität Ihrer Daten beeinträchtigen können.

Überprüfen und Auswechseln des Spülgasfilters

Tauschen Sie den Filter aus, wenn er eine gelbe Farbe angenommen hat, sich die Farbe anderweitig verändert hat oder der Filter mit Fremdstoffen oder Partikeln verunreinigt ist.

HINWEIS

4

Wir empfehlen, dass Sie die Vorrichtungen zur Abdichtung und Trocknung aufrechterhalten und/oder das Gerät stets spülen. Geräteschäden, die entstehen, weil das Gerät nicht gespült wurde oder die Maßnahmen für die Abdichtung und Trocknung nicht aufrechterhalten wurden, sind von der Garantie ausgenommen. Wenn Sie Fragen hierzu haben, setzen Sie sich bitte mit uns in Verbindung.

Abbildung 3-1: Der Spülgasfilter befindet sich in der Kunststoffkappe unterhalb des Manometers.



Hinweis Wenden Sie sich für die Bestellung von Teilen an den technischen Kundendienst.

So wechseln Sie den Spülgasfilter:

- 1. Schließen Sie die Spülgaszufuhr am Hauptventil. Drehen Sie nicht den Durchflussmesser oder Druckregler herunter.
- 2. Entfernen Sie die Kunststoffkappe, in der sich der Filter befindet, und entfernen Sie den Filter. (Sie können beide Teile mit der Hand abschrauben.)



- 3. Setzen Sie einen neuen Filter ein und montieren Sie die Kappe wieder am Manometer.
- 4. Öffnen Sie das Hauptventil, um den Spülgasfluss zum Messgerät zu starten und prüfen Sie auf einen korrekten Spülgasdurchfluss.

[Leerseite]

Kapitel 4

Wartung

Das Nicolet Summit OA FTIR Spektrometer erfordert nur wenig Pflege und keine tägliche Wartung. Durch eine grundlegende Pflege können Sie jedoch eine dauerhafte Spitzenleistung Ihres Spektrometers gewährleisten.

| Wartungsplan für das Summit Spektrometer | |
|---|----|
| Justieren der Spektrometeroptik | 48 |
| Kalibrieren der Laserfrequenz | |
| Reinigen des Spektrometers und Kristalls | 50 |
| Entfernen und Wiedereinsetzen des Kristallaufsatzes | |
| Anbringen der Abdeckung | 54 |
| Auswechseln des Trockenmittels | 59 |
| Austauschen der IR Lichtquelle | 66 |

Wartungsplan für das Summit Spektrometer

Ihr Thermo Scientific Nicolet Summit OA FTIR Spektrometer erfordert nur wenig Pflege und keine tägliche Wartung. Durch eine grundlegende Pflege können Sie jedoch eine dauerhafte Spitzenleistung Ihres Spektrometers gewährleisten. Folgen Sie den Richtlinien in diesem Artikel. Lassen Sie Ihr Spektrometer eingeschaltet, um die bestmögliche Leistung zu erzielen.

HINWEIS

Statische Elektrizität kann die kritischen Bauteile Ihres Spektrometers beschädigen. Beachten Sie die folgenden Empfehlungen, um derartige Beschädigungen zu vermeiden:

- Entladen Sie vor dem Trennen des Netzteils stets die statische Elektrizität, die sich durch Berührungen des Metallgehäuses des Gerätes aufgebaut haben könnte.
- Belassen Sie Ersatz- oder Austauschteile in ihrer Schutzverpackung, bis diese in das Spektrometer eingebaut werden sollen.

Wöchentliche Wartungsmaßnahmen

ÜBERPRÜFEN DER LEISTUNG DES SPEKTROMETERS

Die OMNIC Paradigm-Software umfasst einen Arbeitsablauf zur Leistungsüberprüfung des Spektrometers. Bei dem PV-Arbeitsablauf werden eine Reihe von Standardtests durchgeführt, um die Betriebsbereitschaft des Gerätes zu prüfen und die Genauigkeit Ihrer Daten sicherzustellen.

Wir empfehlen, dass Sie den PV-Arbeitsablauf mindestens einmal wöchentlich ausführen.

REINIGEN DES SUMMIT SPEKTROMETERS

Das Spektrometer darf nur nach den Empfehlungen gereinigt werden. Siehe <u>"Reinigen des Summit Spektrometers"</u>.

Monatliche Wartungsmaßnahmen

ÜBERPRÜFEN DES FEUCHTIGKEITSINDIKATORS

Die optischen Bauteile des Spektrometers können leicht durch übermäßige Luftfeuchtigkeit beschädigt werden. Das Spektrometer ist versiegelt und die enthaltenen optischen Bauteile sind durch zwei Trockenmittelbehälter geschützt, die Feuchtigkeit aufnehmen. Überwachen Sie die Luftfeuchtigkeit des Spektrometers in der OMNIC Paradigm-Software.

HINWEIS

Wir empfehlen, dass Sie das Spektrometer stets geschlossen und mit Trockenmittel versehen halten, oder verwenden Sie eine Spüleinheit, um das Gerät mit Trockenluft oder Stickstoff zu spülen. Geräteschäden, die entstehen, weil das Gerät nicht gespült wurde oder die Maßnahmen für die Abdichtung und Trocknung nicht aufrechterhalten wurden, sind von der Garantie ausgenommen. Wenn Sie Fragen hierzu haben, setzen Sie sich bitte mit uns in Verbindung.

ÜBERPRÜFEN DES SPÜLGASFILTERS

Wenn das Spektrometer mit Stickstoff oder Trockenluft gespült wird, müssen Sie den Spülgasfilter mindestens einmal monatlich überprüfen.

Ersetzen Sie den Filter, wenn er eine gelbe Farbe angenommen hat, sich die Farbe anderweitig verändert hat oder der Filter mit Fremdstoffen oder Partikeln verunreinigt ist. Weitere Informationen finden Sie unter "Überprüfen und Auswechseln des Spülgasfilters".

Abbildung 4-1: Der Spülgasfilter befindet sich in der Kunststoffkappe unterhalb des Manometers.



Justieren der Spektrometeroptik

Die Optik Ihres Spektrometers muss justiert werden, wenn der Arbeitsablauf Nicolet Summit OA - PV-Test fehlgeschlagen ist. Das Justieren des Spektrometers optimiert die Energie, die den Detektor erreicht und maximiert das Detektorsignal. Wenn der Leistungstests fehlschlägt, justieren Sie die Optik des Spektrometers neu, kalibrieren die Laserfrequenz, und führen den Test dann erneut aus. Wenn der Test erneut fehlschlägt, wenden Sie sich an Ihren regionalen technischen Kundendienst.

So justieren Sie die Optik Ihres Spektrometers (Touchscreen-Bedienungsoberfläche):

- 1. Wählen Sie auf der Startseite das Symbol "Diagnose" [1], um die Ansicht "Diagnose" zu öffnen.
- 2. Öffnen Sie die Registerkarte "Justieren" und wählen Sie Justieren.

Das System führt eine automatische Justage des Spektrometers durch und zeigt an, wenn die Justage abgeschlossen ist.

- So justieren Sie die Optik Ihres Spektrometers (Bedienoberfläche):
- 1. Gehen Sie zum Administratorbereich und wählen Sie das Symbol "Diagnose" [100], um die Ansicht "Diagnose" zu öffnen.
- 2. Öffnen Sie die Registerkarte "Justieren" und wählen Sie Justieren.

Das System führt eine automatische Justage des Spektrometers durch und zeigt an, wenn die Justage abgeschlossen ist.

- So justieren Sie die Optik Ihres Spektrometers (Desktop-Bedienungsoberfläche):
- 1. Wählen Daten erfassen > Diagnostik > Spektrometer justieren.
- 2. Klicken Sie im Dialogfeld "Spektrometer justieren" auf Start.

Das System führt eine automatische Justage des Spektrometers durch und zeigt an, wenn die Justage abgeschlossen ist.

Kalibrieren der Laserfrequenz

Die Frequenz des Spektrometers muss kalibriert werden, wenn der Arbeitsablauf Nicolet Summit OA -PV-Test einen fehlgeschlagenen Test anzeigt. Wenn ein Test fehlschlägt, kalibrieren Sie die Laserfrequenz neu und führen den Test dann erneut aus. Wenn der Test erneut fehlschlägt, wenden Sie sich an Ihren regionalen technischen Kundendienst.

So kalibrieren Sie den Laser (Touchscreen-Bedienungsoberfläche):

- 1. Wählen Sie auf der Startseite das Symbol "Diagnose" [1], um die Ansicht "Diagnose" zu öffnen.
- 2. Öffnen Sie die Registerkarte "Laserkalibrierung" und wählen Sie **Kalibrieren**, um den Vorgang zu starten.

Wenn die Kalibrierung des Lasers abgeschlossen hat, zeigt eine Meldung an, ob die Frequenz geändert wurde und bei einer Änderung die entsprechende neue Frequenz.

- So kalibrieren Sie den Laser (Bedienoberfläche):
- 1. Gehen Sie zum Administratorbereich und wählen Sie das Symbol "Diagnose" [100], um die Ansicht "Diagnose" zu öffnen.
- 2. Öffnen Sie die Registerkarte "Laserkalibrierung" und wählen Sie **Kalibrieren**, um den Vorgang zu starten.

Wenn die Kalibrierung des Lasers abgeschlossen hat, zeigt eine Meldung an, ob die Frequenz geändert wurde und bei einer Änderung die entsprechende neue Frequenz.

- So kalibrieren Sie den Laser (Desktop-Bedienungsoberfläche):
- 1. Wählen Sie Daten erfassen > Diagnostik > Laserkalibrierung.
- 2. Klicken Sie in dem Dialogfeld "Laserkalibrierung" auf Start, um den Laser zu kalibrieren.

Wenn das System die Kalibrierung des Lasers abgeschlossen hat, zeigt eine Meldung an, ob die Frequenz geändert wurde und bei einer Änderung die entsprechende neue Frequenz.

Reinigen des Spektrometers und Kristalls

Bevor Sie das Thermo Scientific Nicolet Summit OA Spektrometer reinigen, müssen Sie es abschalten und von der Spannungsversorgung trennen.

Reinigen des Summit Spektrometers

Reinigen Sie die Außenseite des Spektrometers mit einem weichen, sauberen Tuch und einer milden Reinigungslösung wie z. B. Windex. Besprühen Sie das Tuch und wischen Sie ausschließlich die äußere, lackierte Oberfläche ab. Achten Sie darauf, dass keine Reinigungslösung auf die Elektronik hinten am Spektrometer gelangt.

Staub kann sich auf dem Elektronikgehäuse an der Rückseite des Spektrometers ansammeln und potenziell mit der abgestrahlten Wärme reagieren. Das kann sich negativ auf die Lebensdauer der elektronischen Bauteile auswirken.

Blasen Sie Staub, der sich auf der Rückseite des Spektrometers angesammelt hat, mit Druckluft weg. Verwenden Sie keine Flüssigkeit, um den Staub von der Rückseite des Spektrometers zu entfernen.

VORSICHT



Gefahren durch Stromschläge sind zu vermeiden.Lassen Sie keine Flüssigkeiten in das Netzteil oder an die Rückseite des Gerätes gelangen, um Stromschläge zu vermeiden.

HINWEIS

Verwenden Sie keine scharfen Reinigungsmittel, Lösungsmittel, Chemikalien oder Scheuermittel, denn diese können die Oberfläche beschädigen.

Reinigen des ATR-Kristallaufsatzes

Verwenden Sie ein Tuch oder Papiertuch, das sauber, trocken und nicht scheuernd ist, um das ATR-Kristall zu reinigen.

WARNUNG





Zinkselenid (ZnSe) ist giftig. Informationen zur Handhabung und zu Expositionskontrollen entnehmen Sie bitte dem ZnSe-Sicherheitsdatenblatt des Herstellers unter www.specac.com.

Tragen Sie beim Reinigen oder Austauschen des Kristallaufsatzes Schutzhandschuhe.

- 1. Tupfen Sie Flüssigkeiten mit einem sauberen Tuch oder Papiertuch vom Kristall ab. Entsorgen Sie das Papiertuch.
- 2. Tragen Sie mit einer Transferpipette einige Tropfen Heptan auf das Kristall und den umliegenden Bereich auf.
- 3. Wischen Sie das Kristall mit einem neuen, sauberen Papiertuch der Länge nach ab, um alle Flüssigkeiten zu entfernen. Wischen Sie den Bereich um die Kristallmulde ab.



4. Wiederholen Sie den Vorgang bei Bedarf und verwenden Sie abschließend immer ein sauberes, trockenes Tuch oder Papiertuch.

Entfernen und Wiedereinsetzen des Kristallaufsatzes

Der Kristallaufsatz sollte alle 12 bis 18 Monate oder nach Bedarf ausgetauscht werden.

WARNUNG Jas Einatmen und Verschlucken von Giftstoffen ist zu vermeiden. Zinkselenid (ZnSe) ist giftig. Informationen zur Handhabung und zu Expositionskontrollen entnehmen Sie bitte dem ZnSe-Sicherheitsdatenblatt des Herstellers unter www.specac.com. Tragen Sie beim Reinigen oder Austauschen des Kristallaufsatzes Schutzhandschuhe.

HINWEIS

Achten Sie darauf, dass nichts in das Spektrometer fällt, wenn die Abdeckung abgenommen ist.

ERFORDERLICHE MATERIALIEN:

• 3/32-Zoll-Innensechskantschlüssel

Entfernen des Kristallaufsatzes

- 1. Entfernen Sie Staubabdeckung vom Kristall.
- 2. Lösen Sie die drei Schrauben, mit denen der Kristallaufsatz am Spektrometer befestigt ist, mit einem 3/32-Zoll-Innensechskantschlüssel vollständig.



- 3. Heben Sie den Kristallaufsatz in einer geraden Bewegung nach oben vom Spektrometer ab.
- 4. Reinigen Sie den Bereich um den Flansch herum unter der Stelle, an der der Aufsatz befestigt ist.

Wiedereinsetzen des Kristallaufsatzes

- 1. Bringen Sie den neuen Aufsatz über dem Flansch in Position.
- 2. Setzen Sie die Unterlegscheiben auf die Schrauben auf, und setzen Sie die Schrauben in die drei Öffnungen ein. Die Schrauben müssen gerade in die Öffnungen eingesetzt werden.
- 3. Drücken Sie auf den Kristallaufsatz und ziehen Sie eine Schraube auf jeder Seite des Aufsatzes fest. Ziehen Sie dann die dritte Schraube fest, die sich am nächsten zur Rückseite des Spektrometers befindet. Ziehen Sie alle Schrauben fingerfest an.

Wenn sich die Schraube ungehindert drehen lässt, ziehen Sie die Schraube wieder etwas heraus und setzen sie erneut ein, um die Ausrichtung zu korrigieren.

Anbringen der Abdeckung

ERFORDERLICHE WERKZEUGE:

Bevor Sie beginnen, benötigen Sie Folgendes:

- 7/64-Innensechskantschlüssel
- 3/32-Innensechskantschlüssel
- So bringen Sie die Spektrometerabdeckung an:
- 1. Schalten Sie das Spektrometer aus.
 - a. Drücken Sie zum Ausschalten des Spektrometers die Netztaste.
 - b. Ziehen Sie das Netzkabel ab.
 - c. Schalten Sie ggf. die Spülung ab, trennen Sie die Spülleitung vom Spektrometer.
- 2. Entfernen Sie den Kristallaufsatz.
 - a. Entfernen Sie Staubabdeckung vom Kristall.
 - b. Lösen Sie die drei Schrauben vollständig, mit denen der Kristallaufsatz am Spektrometer befestigt ist.



- c. Heben Sie den Kristallaufsatz in einer geraden Bewegung nach oben vom Spektrometer ab.
- d. Reinigen Sie den Bereich um den Flansch herum unter der Stelle, an der der Aufsatz befestigt ist.
- 3. Nehmen Sie die Geräteabdeckung ab.
 - a. Neigen Sie das Spektrometer vorsichtig auf seine Rückseite, sodass es auf den Rippen des Kühlkörpers zum Liegen kommt.
 - b. Lösen Sie mit einem 7/64-Innensechskantschlüssel die vier Schrauben, mit denen die Abdeckung an der Grundplatte des Geräts befestigt ist.



- c. Stellen Sie das Spektrometer vorsichtig wieder auf die Füße. Die vier Schrauben fallen heraus, wenn Sie das Spektrometer wieder auf die Füße stellen. Achten Sie darauf, dass sie nicht verloren gehen.
- d. Heben Sie die Abdeckung in einer geraden Bewegung nach oben ab.



- 4. Bringen Sie die Geräteabdeckung wieder an.
 - a. Bringen Sie die Abdeckung vorsichtig auf der Grundplatte des Geräts in Position.
 - b. Neigen Sie das Spektrometer auf seine Rückseite, sodass es auf den Rippen des Kühlkörpers zum Liegen kommt.
 - c. Setzen Sie die vier Schrauben ein, die Sie zuvor entfernt haben. Ziehen Sie alle Schrauben fest.
 - d. Stellen Sie das Spektrometer wieder auf die Füße.
- 5. Setzen Sie den Kristallaufsatz wieder ein.
 - a. Bringen Sie den Aufsatz über dem Flansch in Position.
 - b. Setzen Sie die Unterlegscheiben auf die Schrauben auf, und setzen Sie die Schrauben in die drei Öffnungen ein. Die Schrauben müssen gerade in die Öffnungen eingesetzt werden.
 - c. Drücken Sie auf den Kristallaufsatz und ziehen Sie eine Schraube auf jeder Seite des Aufsatzes fest. Ziehen Sie dann die dritte Schraube fest, die sich am nächsten zur Rückseite des

Spektrometers befindet. Ziehen Sie alle Schrauben fingerfest an.

Wenn sich die Schraube ungehindert drehen lässt, ziehen Sie die Schraube wieder etwas heraus und setzen sie erneut ein, um die Ausrichtung zu korrigieren.

- 6. Einschalten des Spektrometers
 - a. Schließen Sie das Netzkabel an die Netzsteckdose an und schalten Sie das Gerät am Netzschalter ein.
 - b. Schließen Sie ggf. die Spülleitung wieder am Gerät an, und schalten Sie die Spüleinheit ein.
- 7. Überprüfen der Spektrometerleistung
 - a. Justieren Sie die Spektrometeroptik.
 - i. Wählen Sie im Dashboard der OMNIC Paradigm-Software Daten erfassen > Diagnostik
 > Spektrometer justieren.
 - ii. Wählen Sie **Start**, um mit der Justierung zu beginnen. Bei erfolgreichem Abschluss der Justierung wird auf dem Bildschirm "Justierung abgeschlossen" angezeigt.
 - b. Kalibrieren Sie den Laser.
 - i. Wählen Sie Daten erfassen > Diagnostik > Laserkalibrierung.
 - ii. Wählen Sie Start, um mit der Kalibrierung zu beginnen.

Die Laserkalibrierung nimmt einige Minuten in Anspruch. Nach Abschluss der Kalibrierung zeigt das Display die Meldung "Messung abgeschlossen" an und meldet ggf. Änderungen in der Laserfrequenz.

Das Spektrometer braucht bis zu 12 Stunden, um ein thermisches Gleichgewicht zu erreichen. Um die bestmöglichen Ergebnisse zu erzielen, führen Sie die Ausrichtung und Kalibrierung nach den ersten 12 Betriebsstunden erneut aus.

- c. Führen Sie den Arbeitsablauf Nicolet Summit OA PV-Test aus. Im Rahmen des Arbeitsablaufs der Leistungsüberprüfung wird ein Referenzmaterial verwendet und eine Reihe von Standardtests durchgeführt, um die Betriebsbereitschaft des Spektrometers sowie die Genauigkeit der Messdaten zu überprüfen.
 - i. Kehren Sie zum Dashboard zurück.
 - Führen Sie einen Bildlauf zum Bereich "Arbeitsabläufe" durch und wählen Sie Nicolet Summit OA - PV-Test.

iii. Klicken Sie mit der rechten Maustaste auf den Arbeitsablauf und wählen Sie **Ausführen**, um den Arbeitsablauf zu starten.

Auswechseln des Trockenmittels

Überwachen Sie die Luftfeuchtigkeit des Spektrometers mithilfe der OMNIC Paradigm-Software. Wechseln Sie das Trockenmittel nach Bedarf oder den Richtlinien Ihrer Organisationen aus.

Das Auswechseln des Trockenmittels erfordert das Abnehmen der Geräteabdeckung, das Herausnehmen der Trockenmittelkartusche und das Auswechseln des Trockenmittelbehälters und des O-Rings.

ERFORDERLICHE WERKZEUGE:

Bevor Sie beginnen, benötigen Sie Folgendes:

- 7/64-Innensechskantschlüssel
- Trockenmittel Ersatzkit
- 3/32-Innensechskantschlüssel
- Einen #2 Phillips Kreuzschlitzschraubendreher
- Handschuhe, Fingerlinge oder Labortücher (zur Handhabung des Feuchtigkeitsindikators)

WARNUNG

Explosionsgefahren sind zu vermeiden.



Bevor Sie die Geräteabdeckung entfernen und das Trockenmittelfach öffnen, schalten Sie das Spektrometer aus, ziehen den Netzstecker und entfernen alle Proben aus dem Gerät.

Das Eindringen von entzündlichen Flüssigkeiten und Gasen in das Trockenmittelfach stellt eine Explosionsgefahr dar. Sollte dieser Fall eintreten, setzen Sie sich umgehend mit uns in Verbindung. Das Gerät darf nicht an die Spannungsversorgung angeschlossen werden, bis das Problem behoben wurde.

HINWEIS

Achten Sie darauf, dass keine Gegenstände in das Gerät fallen, während die Trockenmittelabdeckung geöffnet ist.

WARNUNG



Das Trockenmittel darf nur gegen ein von uns geliefertes Ersatzteil ausgetauscht werden.

So wechseln Sie das Trockenmittel aus:

- 1. Schalten Sie das Spektrometer aus.
 - a. Drücken Sie zum Ausschalten des Spektrometers die Netztaste.
 - b. Ziehen Sie das Netzkabel ab.
 - c. Schalten Sie ggf. die Spülung ab, trennen Sie die Spülleitung vom Spektrometer.

Hinweis Sie können entweder eine Trockenmittelkartusche mit oder ohne Spülanschluss verwenden, um die erforderliche Luftfeuchtigkeit aufrechtzuerhalten. Die unten stehenden Abbildungen zeigen eine Trockenmittelkartusche, die nicht gespült wird. Die Anweisungen für den Austausch der Trockenmittelbehälter sind für beide Kartuschenversionen gleich.

- 2. "Entfernen des Kristallaufsatzes"
- 3. Nehmen Sie die Geräteabdeckung ab.
 - a. Neigen Sie das Spektrometer vorsichtig auf seine Rückseite, sodass es auf den Rippen des Kühlkörpers zum Liegen kommt.
 - b. Lösen Sie mit einem 7/64-Innensechskantschlüssel die vier Schrauben, mit denen die Abdeckung an der Grundplatte des Geräts befestigt ist.



- c. Stellen Sie das Spektrometer vorsichtig wieder auf die Füße. Die vier Schrauben fallen heraus, wenn Sie das Spektrometer wieder auf die Füße stellen. Achten Sie darauf, dass sie nicht verloren gehen.
- d. Heben Sie die Abdeckung in einer geraden Bewegung nach oben ab.



- 4. Entfernen der Trockenmittelkartusche
 - a. Lösen Sie die beiden Schrauben mit dem #2 Philipps Kreuzschlitzschraubenzieher von der Trockenmittelkartusche, und ziehen Sie diese gerade nach oben aus dem Spektrometer heraus.



b. Entfernen und entsorgen Sie den großen O-Ring.



Hinweis Achten Sie darauf, dass der O-Ring nicht in die Trockenmittelkammer fällt.

5. Entfernen des Trockenmittelbehälters und ersetzen des Feuchtigkeitsindikators

WARNUNG



Der Inhalt der Trockenmittelbehälter (Aluminiumsilikat Molekularsieb) kann bei Verschlucken schädlich sein. Wenn Sie einen verbrauchten Trockenmittelbehälter entsorgen möchten, achten Sie darauf, dass die Entsorgung den Vorschriften entspricht.

a. Lösen Sie die acht Schrauben, mit denen die zwei gesättigten Trockenmittelbehälter an der Trockenmittelkartusche gehalten werden, mit dem 0,05-Zoll-Innensechskantschlüssel (etwa zwei Umdrehungen). Nehmen Sie die Behälter heraus.



b. Drehen Sie die Trockenmittelkartusche um, und entfernen Sie den alten Feuchtigkeitsindikator vom Fenster. Entsorgen Sie den alten Feuchtigkeitsindikator.

HINWEIS

Verwenden Sie beim Arbeiten mit dem neuen Feuchtigkeitsindikator immer Schutzhandschuhe, Fingerlinge oder Labortücher. Der Indikator kann sich durch Fett oder Feuchtigkeit von der Haut verfärben.

- 6. Einsetzen der neuen Trockenmittelbehälter
 - ä. Öffnen Sie das versiegelte Paket mit dem neuen Trockenmittel (muss versiegelt sein, um die Wirksamkeit des Trockenmittels zu bewahren), und entfernen Sie den neuen Feuchtigkeitsindikator.
 - b. Drehen Sie die Trockenmittelkartusche um und drücken Sie den blauen Indikator auf das Fenster, so dass die flachen Kanten des Indikators mit den Rändern der Kartusche ausgerichtet sind. Der Text sollte durch das Fenster sichtbar sein, wenn Sie die rechte Seite der Kartusche nach oben drehen.

HINWEIS

Acht Sie darauf, dass der Indikator fest gegen das Fenster gedrückt ist, so dass er sich nicht in die Trockenmittelkammer lösen kann, wenn Sie die Kartusche wieder einsetzen.

- c. Setzen Sie die Behälter in den Schlitz der Trockenmittelkartusche ein, und ziehen Sie die acht Schrauben (ungefähr 2 Umdrehungen) fest, bis sie bündig mit dem Trockenmittelkartuschen Schlitz abschließen.
- 7. Einsetzen der Trockenmittelkartusche
 - a. Nehmen Sie den neuen O-Ring aus dem Paket und montieren Sie ihn an der Grundplatte des Messgerätes.

Drücken (besser nicht rollen) Sie den neuen O-Ring in die richtige Position, um sicherzustellen, dass er korrekt in der Nut sitzt.

b. Setzen Sie Trockenmittelkartusche so in das Spektrometer ein, dass der Text auf dem Feuchtigkeitsindikator von einer Person gelesen werden kann, die vor dem Spektrometer steht.

Stellen Sie sicher, dass die Kartusche ordnungsgemäß über dem O-Ring sitzt und ziehen Sie dann die beiden Schrauben mit dem #2 Phillips Kreuzschlitzschraubenzieher fest.

- 8. Bringen Sie die Geräteabdeckung wieder an.
 - a. Bringen Sie die Abdeckung vorsichtig auf der Grundplatte des Geräts in Position.
 - b. Neigen Sie das Spektrometer auf seine Rückseite, sodass es auf den Rippen des Kühlkörpers zum Liegen kommt.
 - c. Setzen Sie die vier Schrauben ein, die Sie zuvor entfernt haben. Ziehen Sie alle Schrauben fest.
 - d. Stellen Sie das Spektrometer wieder auf die Füße.
 - e. "Wiedereinsetzen des Kristallaufsatzes".
- 9. Einschalten des Spektrometers
 - a. Schließen Sie das Netzkabel an die Netzsteckdose an und schalten Sie das Gerät am Netzschalter ein.
 - b. Schließen Sie ggf. die Spülleitung wieder am Gerät an, und schalten Sie die Spüleinheit ein.
- 10. Überprüfen der Spektrometerleistung

- a. Justieren Sie die Spektrometeroptik.
 - Wählen Sie im Dashboard der OMNIC Paradigm-Software Daten erfassen > Diagnostik
 > Spektrometer justieren.
 - ii. Wählen Sie **Start**, um mit der Justierung zu beginnen. Bei erfolgreichem Abschluss der Justierung wird auf dem Bildschirm "Justierung abgeschlossen" angezeigt.
- b. Kalibrieren Sie den Laser.
 - i. Wählen Sie Daten erfassen > Diagnostik > Laserkalibrierung.
 - ii. Wählen Sie Start, um mit der Kalibrierung zu beginnen.

Die Laserkalibrierung nimmt einige Minuten in Anspruch. Nach Abschluss der Kalibrierung zeigt das Display die Meldung "Messung abgeschlossen" an und meldet ggf. Änderungen in der Laserfrequenz.

Das Spektrometer braucht bis zu 12 Stunden, um ein thermisches Gleichgewicht zu erreichen. Um die bestmöglichen Ergebnisse zu erzielen, führen Sie die Ausrichtung und Kalibrierung nach den ersten 12 Betriebsstunden erneut aus.

- c. Führen Sie den Arbeitsablauf Nicolet Summit OA PV-Test aus. Im Rahmen des Arbeitsablaufs der Leistungsüberprüfung wird ein Referenzmaterial verwendet und eine Reihe von Standardtests durchgeführt, um die Betriebsbereitschaft des Spektrometers sowie die Genauigkeit der Messdaten zu überprüfen.
 - i. Kehren Sie zum Dashboard zurück.
 - ii. Führen Sie einen Bildlauf zum Bereich "Arbeitsabläufe" durch und wählen Sie Nicolet Summit OA - PV-Test.
 - iii. Klicken Sie mit der rechten Maustaste auf den Arbeitsablauf und wählen Sie **Ausführen**, um den Arbeitsablauf zu starten.

Austauschen der IR Lichtquelle

Die IR Lichtquelle ist leicht von der Unterseite des Spektrometers aus zugänglich und kann ohne Entfernen der Spektrometerabdeckung ausgetauscht werden.

ERFORDERLICHE WERKZEUGE UND MATERIALIEN:

- Ein #1 Kreuzschlitzschraubenzieher
- IR Lichtquelle Ersatzkit

VORSICHT



Schalten Sie das Spektrometer aus, bevor Sie die IR Lichtquelle austauschen. Die IR Lichtquelle darf nur gegen ein von uns geliefertes Ersatzteil ausgetauscht werden.

So tauschen Sie die IR Lichtquelle aus:

- 1. Ausschalten des Spektrometers und Abziehen des Netzkabels
 - a. Schalten Sie das Spektrometer am Netzschalter aus und ziehen Sie das Netzkabel ab.
 - b. Ziehen Sie alle Kabel vom Spektrometer ab, beispielsweise das Netzwerkkabel und die USB Kabel.
 - c. Wenn das Spektrometer gespült wird, trennen Sie die Spülleitungen vom Spektrometer. Weitere Informationen finden Sie unter <u>"Installieren und Warten einer Spüleinheit"</u>.

VORSICHT



d.

Brandgefahren sind zu vermeiden.

Die IR Lichtquelle wird beim Arbeiten mit dem Gerät sehr heiß. Warten Sie mindestens 10 Minuten nach dem Abschalten des Gerätes, bevor Sie mit dem nächsten Schritt fortfahren.

Thermo Scientific

- 2. Ausbau der IR Lichtquelle
 - a. Lösen Sie die drei Schrauben, mit denen die IR Lichtquelle fixiert ist, mithilfe eines #1-Kreuzschlitzschraubenzieher.



b. Schwenken Sie den Griff an der Lichtquelle nach oben. Halten Sie das Spektrometer mit einer Hand und ziehen Sie fest an dem Griff, um die IR Lichtquelle gerade aus dem Spektrometer herauszuziehen.



HINWEIS

Die Komponenten der Lichtquelle dürfen niemals mit den bloßen Fingern berührt werden. Hautfett und andere Ablagerungen verkürzen die Lebensdauer der Lichtquelle. Verwenden Sie bei der Handhabung einer IR Lichtquelle immer saubere Fingerlinge, Handschuhe oder saubere Labortücher.

- 3. Einsetzen der neuen IR Lichtquelle
 - a. Achten Sie darauf, dass die drei Schrauben gerade sind und schieben Sie die IR Lichtquelle in die Aussparung, bis die Halterung der IR-Quelle bündig mit dem Spektrometer abschließt.
 - b. Ziehen Sie die Schrauben fest, um die IR Lichtquelle zu fixieren, und schwenken Sie den Griff zurück in die Parkposition.
- 4. Anschließen und Einschalten des Spektrometers

- a. Stellen Sie das Messgerät vorsichtig wieder in die aufrechte Position und schließen Sie alle zuvor entfernten Kabel wieder an.
- b. Wechseln Sie das Trockenmittel aus. Anweisungen hierzu finden Sie unter <u>"Auswechseln des</u> <u>Trockenmittels"</u>.

Da die internen Bauteile der Optik beim Austauschen der IR Lichtquelle der Atmosphäre ausgesetzt werden, sollte nach dem Austausch der IR Lichtquelle auch jedes Mal das Trockenmittel gewechselt werden.

WARNUNG



Gefahren durch Stromschläge sind zu vermeiden.

Um eine gute Erdverbindung zu gewährleisten und die Gefahr von Stromschlägen zu vermeiden, verwenden Sie keine Steckdose, die mit einer Erdungsleitung verbunden ist. Die Erdung muss über ein nicht stromführendes Kabel am Hauptverteilerkasten erfolgen.

- c. Schließen Sie das Netzkabel an die Netzsteckdose an und schalten Sie das Gerät am Netzschalter ein.
- 5. Überprüfen der Spektrometerleistung
 - a. Justieren Sie die Spektrometeroptik.
 - i. Wählen Sie im Dashboard der OMNIC Paradigm-Software **Daten erfassen > Diagnostik > Spektrometer justieren**.
 - ii. Wählen Sie **Start**, um mit der Justierung zu beginnen. Bei erfolgreichem Abschluss der Justierung wird auf dem Bildschirm "Justierung abgeschlossen" angezeigt.
 - b. Kalibrieren Sie den Laser.
 - i. Wählen Sie Daten erfassen > Diagnostik > Laserkalibrierung.
 - ii. Wählen Sie Start, um mit der Kalibrierung zu beginnen.

Die Laserkalibrierung nimmt einige Minuten in Anspruch. Nach Abschluss der Kalibrierung zeigt das Display die Meldung "Messung abgeschlossen" an und meldet ggf. Änderungen in der Laserfrequenz.

Das Spektrometer braucht bis zu 12 Stunden, um ein thermisches Gleichgewicht zu erreichen. Um die bestmöglichen Ergebnisse zu erzielen, führen Sie die Ausrichtung und Kalibrierung nach den ersten 12 Betriebsstunden erneut aus.

- c. Führen Sie den Arbeitsablauf Nicolet Summit OA PV-Test aus. Im Rahmen des Arbeitsablaufs der Leistungsüberprüfung wird ein Referenzmaterial verwendet und eine Reihe von Standardtests durchgeführt, um die Betriebsbereitschaft des Spektrometers sowie die Genauigkeit der Messdaten zu überprüfen.
 - i. Kehren Sie zum Dashboard zurück.
 - ii. Führen Sie einen Bildlauf zum Bereich "Arbeitsabläufe" durch und wählen Sie Nicolet Summit OA - PV-Test.
 - iii. Klicken Sie mit der rechten Maustaste auf den Arbeitsablauf und wählen Sie **Ausführen**, um den Arbeitsablauf zu starten.
[Leerseite]