Nicolet Summit OA

Spettrometro FTIR



GUIDA PER L'UTENTE

269-346500 Revisione B Aprile 2021

© 2020 Thermo Fisher Scientific Inc. Tutti i diritti riservati.

Per il Supporto tecnico, contattare il sito: www.thermofisher.com

Al momento dell'acquisto, Thermo Fisher Scientific Inc. fornisce ai propri clienti il presente documento che dovrà essere utilizzato durante il funzionamento del prodotto. Il presente documento è tutelato dai diritti d'autore e la sua riproduzione, intera o parziale, è severamente vietata se non autorizzata per iscritto da Thermo Fisher Scientific Inc.

Il contenuto del presente documento è soggetto a modifiche senza preavviso. Tutte le informazioni tecniche contenute nel presente documento hanno puro scopo di riferimento. Le configurazioni e le specifiche del sistema riportate nel presente documento sostituiscono e annullano tutte le precedenti informazioni ricevute dall'acquirente.

Thermo Fisher Scientific Inc. non garantisce che il presente documento sia completo, accurato o privo di errori e non si assume alcuna responsabilità per errori, omissioni, danni o perdite che possano derivare dall'uso del documento, anche seguendo correttamente le istruzioni che vi sono contenute.

Il presente documento non rientra in un contratto di vendita tra Thermo Fisher Scientific Inc. e l'acquirente. Il documento non regola né modifica i Termini e le condizioni di vendita; in caso di conflitto tra le informazioni contenute nei due documenti, si dovrà fare riferimento ai Termini e alle condizioni di vendita.

Solo per ricerca. Questo strumento o accessorio non è un dispositivo medico e non è destinato all'uso a scopo di prevenzione, diagnosi, trattamento o cura di malattie.



Pericolo di esplosione o incendio.

Questo strumento o accessorio non è destinato all'uso in atmosfera esplosiva.

Sommario

Benvenuto	1
Convenzioni utilizzate	2
Ordine delle parti	3
Contattare la nostra azienda	3

Iniziare con lo spettrometro Nicolet Summit OA	5
Prima della consegna dello spettrometro Summit	6
Sicurezza personale con lo spettrometro Summit	9
Disimballaggio dello spettrometro e operazioni preliminari	
Backup e ripristino dei dati	
Installazione di un adattatore USB wireless	
Informazioni sulla barra luminosa dello spettrometro	
Visualizzazione dei dati sulla app OMNIC Anywhere	31

Opzioni e accessori	33
Installazione e manutenzione di un kit di spurgo	34

Manutenzione	45
Pianificazione della manutenzione per lo spettrometro Summit	46
Allineamento delle ottiche dello spettrometro	48
Calibrazione della frequenza del laser	49
Pulizia dello strumento e del cristallo	50
Rimozione e riposizionamento della piastra del cristallo	52
Riposizionamento del coperchio	54
Sostituzione dell'essiccante	
Sostituzione della sorgente	64

Capitolo 1

Benvenuto

Lo spettrometro infrarosso a trasformata di Fourier (FTIR) Nicolet Summit OA di Thermo Scientific permette di eseguire analisi chimiche di campioni acquisendo i dati nel range spettrale del medio infrarosso. Il sistema dispone di funzioni di verifica integrate, di una potente suite software e di numerose altre funzioni che consentono di semplificare l'acquisizione dei dati. È possibile installare autonomamente l'hardware opzionale ed eseguire senza richiedere interventi diverse procedure di assistenza e manutenzione. Il presente documento, o altra documentazione fornita, contiene tutte le informazioni necessarie.

Avere cura di leggere la guida alla sicurezza fornita in dotazione con il sistema prima di utilizzare lo strumento.

Nota È consigliabile mantenere sempre isolato ed essiccato e/o spurgare lo strumento. I danni all'apparecchiatura causati dal mancato isolamento ed essiccazione e/o spurgo non sono coperti dalla garanzia. In caso di domande in merito, rivolgersi alla nostra azienda.

Convenzioni utilizzate

Precauzioni di sicurezza e altre informazioni importanti presentate nel seguente formato:

PERICOLO



Pericolo.Indica una situazione potenzialmente pericolosa che, se non evitata, causerà morte o lesioni gravi.

AVVERTENZA



Pericolo. Indica una situazione potenzialmente pericolosa che, se non evitata, potrebbe provocare la morte o lesioni gravi.

ATTENZIONE



Pericolo. Indica una situazione pericolosa che, se non evitata, potrebbe provocare lesioni di entità lieve o moderata.

AVVISO

Attenersi alle istruzioni evidenziate con questo simbolo per evitare di danneggiare l'hardware del sistema o di perdere dati.

Nota Contiene informazioni aggiuntive utili.

Benvenuto

Ordine delle parti

Per ordinare parti, vi preghiamo di contattare la nostra azienda nei Contatti.

Per inviarci lo strumento o un accessorio per la riparazione, potete prima contattarci telefonicamente o via e-mail per qualsiasi esigenza ai fini della spedizione o per richiedere altre istruzioni.

Contattare la nostra azienda

Per il Supporto tecnico, contattare il sito: www.thermofisher.com

[Pagina lasciata intenzionalmente vuota]

Iniziare con lo spettrometro Nicolet Summit OA

Per iniziare a utilizzare il vostro nuovo spettrometro Nicolet Summit OA, esaminare i requisiti di preparazione del sito e di sicurezza prima che lo strumento venga consegnato, disimballato e configurato e che vi vengano collegati ulteriori accessori.

Prima della consegna dello spettrometro Summit	6
Sicurezza personale con lo spettrometro Summit	
Disimballaggio dello spettrometro e operazioni preliminari	15
Backup e ripristino dei dati	
Installazione di un adattatore USB wireless	
Informazioni sulla barra luminosa dello spettrometro	
Visualizzazione dei dati sulla app OMNIC Anywhere	

Prima della consegna dello spettrometro Summit

Preparare un'area di lavoro adatta prima della consegna dello spettrometro per garantire i dati più accurati e le prestazioni più affidabili a lungo termine.

Questa sezione spiega cosa fare all'arrivo dello spettrometro ed introduce i fattori ambientali ed elettrici che possono influire sulle prestazioni dello spettrometro. Per una spiegazione più esauriente di tali requisiti, consultare le informazioni sul sito e la sicurezza per lo spettrometro.

Disimballaggio dello spettrometro

All'arrivo dello strumento, prima di aprire la scatola, eseguire due importanti passaggi:

• Controllare l'esterno dell'imballo di spedizione per verificare che non abbia subito danni

Se si notano segni di danni, potete contattarci o rivolgervi alla filiale Thermo Fisher Scientific locale per ricevere istruzioni.

• Attendere che lo spettrometro raggiunga la temperatura ambiente

All'interno della scatola, lo spettrometro è sigillato in un sacco di plastica in modo che si mantenga asciutto. **Prima di aprire il sacco, attendere 24 ore che lo spettrometro raggiunga la temperatura ambiente.** Aprendo il sacco prima che lo spettrometro si sia stabilizzato, si potrebbe formare una condensa dannosa per le ottiche interne e causa di danni permanenti.

La garanzia non copre danni dovuti a movimentazioni non appropriate o alla rimozione del sacco di plastica sigillato prima che lo strumento abbia raggiunto la temperatura ambiente.

Preparazione dell'area di lavoro

Prima della consegna dello strumento, verificare che l'area di lavoro sia in grado di sostenere adeguatamente il peso dello spettrometro. Oltre a uno spazio sufficiente per lo spettrometro, prendere in considerazione i diversi requisiti ambientali ed elettrici richiesti.

Dimensioni dello spettrometro

Pur avendo lo spettrometro Summit un ingombro piuttosto ridotto, assicurarsi di lasciare attorno allo strumento uno spazio sufficiente per la dissipazione del calore dagli sfiati e un agevole accesso alle porte, all'alimentazione e ai cavi.

- Peso dello strumento Summit: 12,6 Kg
- Dimensioni (L x A x P): 34 cm x 24 cm x 32 cm

Fattori ambientali

Lo spettrometro Summit è uno strumento robusto progettato per l'uso in numerosi ambienti. Per ottenere le prestazioni migliori è tuttavia opportuno tenerlo in un ambiente relativamente privo di polvere e con un tasso di umidità basso. Lo spettrometro garantisce un funzionamento affidabile a temperature fra 15 e 35 °C ma per ottenere prestazioni ottimali è consigliabile tenerlo a una temperatura di 20-22 °C.

L'umidità può causare formazione di condensa al suo interno e danneggiare i componenti interni. Alcune precauzioni possono generalmente consentire di proteggere lo strumento dall'umidità:

- Lasciare l'essiccante all'interno dello strumento anche quando viene riposto per la conservazione.
- Evitare repentine variazioni di temperatura.
 - Tenere lo strumento lontano da fonti di aria fredda o calda, come ad esempio in prossimità degli sfiati di impianti di riscaldamento e condizionamento oppure vicino a grandi finestre.

Se viene tenuto in un ambiente particolarmente umido, prendere in considerazione di installare un kit di gas di spurgo.

Requisiti elettrici

L'alimentazione dello spettrometro deve essere fornita da fonti di alimentazione ininterrotta dedicate e prive dei seguenti disturbi:

- Cadute di tensione
- Picchi transitori
- Variazioni di frequenza
- Altri disturbi di rete

Se si sospettano problemi di alimentazione si consiglia di eseguire un controllo della qualità della tensione. Contattare la nostra azienda o la società elettrica locale per maggiori informazioni.

Specifiche dell'alimentazione elettrica

La tabella seguente riporta le specifiche per l'alimentazione elettrica. Per qualsiasi chiarimento sui requisiti, rivolgersi al responsabile del nostro servizio di assistenza locale.

Iniziare con lo spettrometro Nicolet Summit OA

Requisiti	Specifica
Corrente in ingresso	1,5 A max
Tensione in ingresso	Da 100 a 240 V c.a.
Frequenza di linea	da 50 a 60
Disturbi sulla linea	Abbassamenti di tensione, sovracorrenti momentanee e altri disturbi sulla rete non devono superare il 10% della tensione in ingresso
Rumore	meno di 2 V (modo comune) meno di 20 V (modo normale)

Sicurezza personale con lo spettrometro Summit

Pur essendo progettato per essere uno strumento sicuro, prendendo alcune precauzioni è possibile proteggersi da potenziali pericoli che possono insorgere durante il normale utilizzo e manutenzione dello spettrometro Nicolet Summit OA.

ATTENZIONE



Questa guida costituisce un'introduzione ai potenziali pericoli di cui tenere conto ma non è una guida esaustiva. Prima di utilizzare lo strumento consultare le informazioni sul sito e la sicurezza per una descrizione completa di questi potenziali pericoli.

Potenziali pericoli durante il normale utilizzo

Durante il normale utilizzo, la maggior parte dei pericoli deriva dalle seguenti fonti:

- Campioni e solventi potenzialmente pericolosi
- Esposizione al seleniuro di zinco (ZnSe)
- Parti dello spettrometro calde o che emanano calore
- Il laser dello strumento

Comprendendo i potenziali pericoli e prendendo alcune precauzioni è possibile evitare di farsi male o di danneggiare lo strumento.

Campioni e solventi pericolosi

Prendere le opportune precauzioni quando si utilizzano o si programma di misurare campioni o solventi potenzialmente pericolosi quali gas pressurizzati o solventi corrosivi o infiammabili.

VENTILAZIONE APPROPRIATA

Non sono previsti particolari requisiti per quanto riguarda la ventilazione dello spettrometro ma durante alcuni tipi di analisi potrebbe essere necessario impiegare una ventilazione aggiuntiva. Assicurare una ventilazione adeguata durante l'analisi di campioni altamente tossici, la dissoluzione dei campioni in solventi che possono interagire con la sorgente a infrarossi o il campionamento di gas infiammabili.

Iniziare con lo spettrometro Nicolet Summit OA

La pirolisi di solventi che contengono idrocarburi alogenati potrebbe produrre acido cloridrico (HCI), acido fluoridrico (HF) o fosgene (COCI₂).

AVVERTENZA



Pericolo di inalazione di sostanze tossiche. L'acido cloridrico, l'acido fluoridrico e il fosgene sono altamente tossici. Se si utilizzano solventi che contengono idrocarburi alogenati, accertarsi che l'area di lavoro sia correttamente ventilata.

SOLVENTI VOLATILI E INFIAMMABILI

La sorgente infrarossa contenuta all'interno dello spettrometro può incendiare campioni e solventi infiammabili e volatili. Adottare le seguenti misure quando si lavora con campioni e solventi infiammabili:

- Lavorare con le finestre del vano campioni installate.
- Assicurare che l'area di lavoro sia opportunamente ventilata con un sistema di ventilazione attivo privo di scintille o altre fonti di accensione e atto ad impedire la raccolta di vapori infiammabili nell'atmosfera circostante lo strumento.
- Non lasciare solventi o campioni infiammabili accanto allo strumento.
- Non lasciare solventi o campioni infiammabili accanto allo strumento nel vano campioni più a lungo del necessario.
- Utilizzare solo aria secca e pulita o azoto per spurgare lo spettrometro.

SOLVENTI CORROSIVI

L'uso di solventi che producono vapori di HCI o HF nel vano campioni può danneggiare gravemente il sistema. Se si utilizzano solventi alogenati, spurgare lo strumento con aria secca e pulita o azoto.

AVVISO

La garanzia non copre i danni all'apparecchiatura causati dal mancato spurgo dello strumento.

MATERIALI A RISCHIO BIOLOGICO O RADIOATTIVI E AGENTI INFETTIVI

Campioni biologici come tessuti, fluidi corporei, agenti infettivi e sangue possono potenzialmente trasmettere malattie infettive. Attenersi ai protocolli del Programma di sicurezza biologica implementati dell'azienda per operare con materiali potenzialmente infettivi.

Materiali tossici

Lo spettrometro Nicolet Summit OA include un cristallo ATR in seleniuro di zinco (ZnSe) e altri strumenti Summit possono utilizzare le finestre del vano campioni in ZnSe opzionali.

AVVERTENZA



Pericolo di inalazione e ingestione di sostanze tossiche.

Il seleniuro di zinco (ZnSe) è tossico. Fare riferimento alla Scheda di dati di sicurezza ZnSe del produttore all'indirizzo www.specac.com per la manipolazione e i controlli dell'esposizione.

Fonti di calore

Alcune parti dello spettrometro possono diventare molto calde durante il normale utilizzo. Prestare attenzione alla sorgente a infrarossi e agli sfiati dello spettrometro.

La superficie rivolta verso l'esterno della sorgente infrarossa si trova sotto lo strumento e può diventare molto calda. Non toccare la parte inferiore dello strumento mentre è in uso o poco dopo l'utilizzo.

Lo sfiato dello spettrometro si trova sul lato sinistro dello strumento. Durante il normale utilizzo, lo spettrometro dissipa aria calda dai componenti interni all'area all'esterno dello spettrometro attraverso questi sfiati. Lasciare spazio sufficiente attorno agli sfiati dello strumento per dissipare l'aria calda.

Sicurezza del laser e delle ottiche

Durante il normale utilizzo dello spettrometro non si è mai esposti a livelli di radiazioni laser non sicuri. In caso di rimozione del coperchio durante una procedura di assistenza si consiglia di prendere particolari precauzioni come l'uso di protezioni oculari. Il tecnico dell'assistenza provvederà a notificare tale necessità.

AVVERTENZA



Pericolo di lesioni personali.

Non fissare mai direttamente il raggio laser o la luce da esso riflessa. Non manomettere mai il laser nemmeno per sostituire un laser difettoso; ciò potrebbe provocare un'esposizione diretta alla luce del laser o all'alta tensione.

Potenziali pericoli durante la manutenzione

Durante gli interventi di manutenzione sullo strumento si potrebbe essere esposti a pericoli diversi rispetto a quelli ai quali si va incontro durante il normale utilizzo. I principali pericoli che possono verificarsi quando si eseguono interventi di manutenzione comprendono lo spurgo dello strumento e l'utilizzo dei suoi componenti interni.

Spurgo dello strumento

In ambienti estremamente umidi, è consigliabile installare una fonte di aria pulita e secca o azoto per spurgare lo spettrometro. Lo spurgo dello strumento può contribuire a proteggere le ottiche interne dal danno causato da un ambiente umido o solventi corrosivi e può garantire risultati più accurati.

Consultare <u>"Installazione e manutenzione di un kit di spurgo"</u> per dettagli sull'acquisto e l'installazione di un kit di spurgo per lo spettrometro.

PERICOLO

Pericolo di incendio ed esplosione.



- Utilizzare solo aria secca o azoto per spurgare lo spettrometro.
- Non utilizzare mai un gas infiammabile, combustibile o tossico per spurgare lo strumento. Il gas di spurgo deve essere privo di olio o altri materiali reattivi. Il calore emesso dalla sorgente o dall'assorbimento del laser potrebbe fare incendiare gas infiammabili o altri materiali reattivi contenuti nel gas di spurgo.

Utilizzo dei componenti interni

Di norma non dovrebbe esserci motivo di rimuovere il coperchio dello strumento o di tentare di sostituire i componenti interni. Tuttavia, se occorre rimuovere il coperchio per procedere alla manutenzione, tenere conto dei rischi di scossa elettrica, ustione o esposizione alla luce laser.

ATTENZIONE



Pericolo di scosse elettriche.

Anche dopo che lo strumento è stato scollegato da tutte le sorgenti di alimentazione elettrica, i condensatori possono mantenere la carica per altri 30 secondi e causare scosse elettriche.

ATTENZIONE



Pericolo di ustioni.

I componenti interni, soprattutto la sorgente infrarossa, possono diventare estremamente caldi durante il normale utilizzo. Spegnere lo strumento e attendere almeno 10 minuti prima di sostituire un componente.

AVVERTENZA



Pericolo di lesioni personali.

- Non fissare mai direttamente il raggio laser o la luce da esso riflessa. Non manomettere mai il laser. Si verrebbe esposti alla radiazione del laser o all'alta tensione.
- Se si eseguono regolazioni del laser o procedure non descritte nei manuali e nelle guide per l'utente, si potrebbe essere esposti a radiazioni pericolose.

SOSTITUZIONE DELL'ESSICCANTE

Nell'aprire il vano dell'essiccante, evitare la penetrazione al suo interno di liquidi o gas infiammabili. Per istruzioni su come sostituire l'essiccante, vedere <u>"Sostituzione dell'essiccante"</u>.

PERICOLO



Pericolo di esplosione.

Prima di aprire il vano dell'essiccante, spegnere lo strumento, scollegare il cavo di alimentazione e rimuovere i campioni dal sistema. La penetrazione di liquidi o gas infiammabili nel vano dell'essiccante potrebbe provocare un'esplosione. In tale evenienza, contattare immediatamente la nostra azienda e non alimentare elettricamente lo strumento finché la condizione non sia stata risolta.

SOSTITUZIONE DELLA SORGENTE IR

La sorgente IR diventa estremamente calda durante l'uso. Se occorre sostituirla, evitare pericoli di ustione ed esplosione.

ATTENZIONE



Pericolo di ustioni.

La sorgente diventa estremamente calda durante il normale utilizzo. Dopo lo spegnimento dello spettrometro, lasciarla sempre raffreddare per almeno 10 minuti prima di cominciare a utilizzarla.

PERICOLO





Prima di togliere la sorgente dallo spettrometro, spegnere lo strumento, scollegare il cavo di alimentazione, staccare le linee di spurgo e rimuovere i campioni dal sistema. La penetrazione di liquidi o gas infiammabili nel vano della sorgente potrebbe provocare un'esplosione. In tale evenienza, contattare immediatamente la nostra azienda e non alimentare elettricamente lo strumento finché la condizione non sia stata risolta.

Riepilogo

Lo spettrometro Nicolet Summit OA è uno strumento sicuro e robusto ma durante l'uso e gli interventi di manutenzione si potrebbe essere esposti a pericoli. Durante il normale utilizzo, fare attenzione nel manipolare campioni e solventi potenzialmente pericolosi ed evitare componenti dello spettrometro caldi o che emettono calore. Durante gli interventi manutentivi, prendere opportune precauzioni per evitare danni o lesioni che potrebbero verificarsi durante lo spurgo dello strumento, la manipolazione dei componenti interni o la sostituzione dell'essiccante.

Disimballaggio dello spettrometro e operazioni preliminari

Il disimballaggio e la configurazione dello spettrometro Nicolet Summit OA di Thermo Scientific richiedono il suo posizionamento nell'area di lavoro, il collegamento e l'accensione nonché l'esecuzione di un breve test di valutazione delle prestazioni.

Prima di cominciare

AVVISO

Attendere 24 ore prima di estrarre lo spettrometro dal sacco di plastica per evitare che la formazione di condensa possa danneggiare le ottiche interne. La garanzia non copre i danni dovuti all'apertura del sacco di plastica sigillato prima che lo strumento abbia raggiunto la temperatura ambiente.

Promemoria: Esaminare le Informazioni sul sito e la sicurezza per assicurarsi che l'area di lavoro sia sicura e idonea per lo strumento.

Disimballaggio e configurazione

Le fasi necessarie per il disimballaggio e la configurazione dello spettrometro variano in relazione alle opzioni acquistate.

1. Disimballare la scatola di spedizione

Estrarre lo spettrometro dalla scatola di spedizione e dal sacco e togliere i restanti materiali di imballaggio.

Lo spettrometro è fornito dotato di un coperchio parapolvere rimovibile sopra il cristallo ATR. Utilizzare il coperchio parapolvere quando si ripone si sposta lo strumento e rimuoverlo durante il normale utilizzo.

- 2. Collegare l'alimentatore allo spettrometro e a una sorgente di alimentazione CA correttamente collegata a terra.
- 3. Collegare il monitor, la tastiera e il mouse.
 - a. Collegare il monitor al connettore Mini DisplayPort dello strumento.
 - b. Collegare la tastiera e il mouse a una delle porte USB dello spettrometro.

Nota Alcuni dispositivi USB e Mini DisplayPort sono noti per interferire con il segnale emesso da dispositivi USB wireless, come le tastiere e i mouse wireless. Se si riscontrano problemi di prestazioni con la tastiera e il mouse wireless, come ritardi o input tardivi o assenti, utilizzare la prolunga USB per posizionare il ricevitore USB wireless più lontano dalle porte USB.

4. Premere l'interruttore di alimentazione dello strumento per accendere lo spettrometro. La barra luminosa dello spettrometro visualizza una luce verde ciclica mentre è in corso la fase di riscaldamento dello strumento e una barra verde fissa quando è pronto per l'uso.

Quando si accende e si spegne lo strumento è possibile percepire il rumore del riferimento interno di polistirene che si sposta. Questo è normale.



Alla prima accensione dello spettrometro, è richiesto l'inserimento di una nuova password. Si consiglia di non immettere una password in questa fase. Procedere senza immettere la password quindi, dopo aver completato la configurazione, gestire gli account e le password sul dispositivo secondo i criteri applicati dalla propria organizzazione.

Nota Se quando si accende il sistema lo schermo è nero, controllare l'alimentazione del monitor, il connettore DisplayPort e l'input del monitor. Consultare la documentazione fornita con il monitor per conoscere i dettagli per la modifica delle impostazioni di input.

5. Aprire il software OMNIC Paradigm.

Il software può impiegare diversi minuti per avviarsi e collegarsi allo strumento. Il software visualizza lo stato di connessione nell'angolo superiore destro della schermata. Quando lo stato evidenzia che si è connessi, procedere al passaggio successivo.



Nota Se si apre un menu nel software prima che lo strumento sia connesso, la funzione Allinea e Calibra sarà disabilitata. Per abilitarla, quando lo strumento è connesso, fare clic sul Dashboard nella barra degli strumenti.

- 6. Verificare le prestazioni dello spettrometro.
 - a. Allineamento delle ottiche dello spettrometro.
 - i. Dall'interfaccia Dashboard andare a Acquisisci dati > Diagnostica > Allinea spettrometro.
 - ii. Selezionare **Avvio** per iniziare l'allineamento. La schermata visualizza "Allineamento completato" quando l'allineamento è terminato con successo.
 - b. Calibrare il laser.
 - i. Andare a Acquisisci dati > Diagnostica > Calibrazione laser.
 - ii. Selezionare Avvio per iniziare la calibrazione.
 - iii. Il completamento della calibrazione laser può richiedere diversi minuti. Terminata la calibrazione, il display visualizza "Misura completata" e indica le modifiche della frequenza laser.

Lo strumento impiega 12 ore per raggiungere l'equilibrio termico. Per conseguire i migliori risultati, rieseguire l'allineamento e la calibrazione dopo le prime 12 ore di funzionamento.

c. Eseguire il workflow di verifica delle prestazioni Nicolet Summit OA .

Il workflow di verifica delle prestazioni utilizza un materiale di riferimento per eseguire una serie di prove standard per verificare il funzionamento dello strumento e garantire l'accuratezza dei dati.

- i. Dal Dashboard dell'interfaccia Desktop, scorrere verso il riquadro Workflow.
- ii. Nel riquadro Workflow, fare clic con il pulsante destro del mouse sul workflow Verifica delle prestazioni (PV) Nicolet Summit OA e selezionare Esegui.

Iniziare con lo spettrometro Nicolet Summit OA



- iii. Seguire le istruzioni visualizzate sullo schermo per completare il workflow.
- iv. Quando viene visualizzata la tabella dei risultati PV, selezionare Fine per completare il workflow e tornare al dashboard o selezionare Misura spettro Squalene per procedere alla porzione dello squalene opzionale del workflow. La misurazione di uno spettro dello squalene costituisce una verifica di riferimento aggiuntiva delle prestazioni del sistema.



Fasi successive

Congratulazioni! Il vostro strumento è pronto per l'utilizzo. Prima di avviare l'acquisizione dei dati occorre tuttavia eseguire il backup del sistema e creare un'unità di ripristino per proteggere il sistema da potenziali problemi futuri. Windows 10 include diversi strumenti integrati di backup e ripristino dei dati.

Per istruzioni per la creazione di un'unità di ripristino e dell'immagine di sistema e per il ripristino del sistema, vedere <u>"Backup e ripristino dei dati"</u>.

Per maggiori informazioni sull'identificazione e l'analisi dei campioni con lo spettrometro, vedere le guide e le esercitazioni di OMNIC Paradigm alla pagina thermofisher.com/ftir-help.

Iniziare con lo spettrometro Nicolet Summit OA

Per commutare all'interfaccia Operatore, andare a Vista / Visualizza > Operatore.

Per istruzioni dettagliate sulla creazione e l'esecuzione di soluzioni personalizzate utilizzando l'editor pacchetti e l'interfaccia Operatore, vedere "Soluzioni personalizzate" nella guida per l'utente del software OMNIC Paradigm.

Backup e ripristino dei dati

Prima di iniziare la raccolta dei dati, prendere alcune precauzioni per proteggere il sistema da eventuali problemi in futuro. Windows 10 include diversi strumenti integrati di backup e ripristino dei dati.

IMPORTANTE! Per garantire la possibilità di ripristinare i dati nel caso in cui il disco rigido o il sistema operativo vengano danneggiati, creare dei supporti di ripristino, creare un'immagine di sistema e stabilire procedure per il backup dei dati.

Creazione di un'unità di ripristino

Creare un'unità di ripristino per potere accedere ad opzioni di avvio avanzate di Windows 10 quali Ripristino immagine di sistema, Ripristino del sistema e Impostazioni di avvio, anche se nell'installazione di Windows vengono danneggiate.

Materiali richiesti

• Unità flash formattata con FAT32 con almeno 8 GB di spazio di archiviazione

AVVISO

Durante questa procedura tutti i dati contenuti nell'unità verranno eliminati. Prima di procedere, accertarsi di avere eseguito il backup di tutti i file che si desidera conservare.

Per creare un'unità di ripristino

- 1. Inserire l'unità flash in una delle porte USB presenti sul lato dello spettrometro.
- 2. Nella casella di ricerca della barra delle applicazioni di Windows 10, cercare "Crea unità di ripristino" quindi selezionare **Crea unità di ripristino**.



3. Alla richiesta di consentire di apportare modifiche al dispositivo, selezionare Sì.

Iniziare con lo spettrometro Nicolet Summit OA

- 4. Assicurarsi di deselezionare la casella di controllo "Esegui il backup dei file di sistema nell'unità di ripristino" e fare clic sul pulsante **Avanti**.
- 5. Selezionare l'unità dall'elenco di quelle disponibili e fare clic sul pulsante Avanti.



- 6. Fare clic su Crea. Al termine del ripristino del sistema, fare clic su Fine per chiudere la finestra.
- 7. Rimuovere l'unità flash in modo sicuro e conservarla in un luogo sicuro. È consigliabile etichettarla come Unità di ripristino Summit.

Per maggiori informazioni sulla creazione di supporti di ripristino, vedere la guida in linea di Microsoft.

Creazione di un'immagine di sistema

Un'immagine di sistema è una foto istantanea dell'intero sistema, comprendente i dati e i file, le applicazioni installate e altre applicazioni e impostazioni. In caso di disco rigido completo o di errore del sistema operativo, utilizzare l'immagine di sistema per ripristinare il dispositivo allo stato in cui era al momento della creazione dell'immagine. La frequenza con cui si desidera aggiornare l'immagine di sistema deve essere valutata soggettivamente ma è consigliabile conservare un'immagine di un'installazione pulita della prima configurazione dello strumento.

A seconda della quantità dei dati copiati nell'immagine di sistema, la procedura potrebbe richiedere tempo. Se viene copiata una grande quantità di dati, il completamento del processo potrebbe richiedere più di un'ora.

Materiali richiesti

SSF formattato da NTFS

Iniziare con lo spettrometro Nicolet Summit OA

Per creare un'immagine di sistema

- 1. Collegare I'SSD a una delle porte USB presenti sullo spettrometro.
- 2. In Windows 10, aprire il Pannello di controllo e selezionare Backup e ripristino (Windows 7).
- 3. Nel pannello a sinistra, selezionare Crea immagine di sistema.

🛞 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. Selezionare Su disco rigido quindi selezionare l'SSD dall'elenco. Fare clic sul pulsante Avanti.

🏥 Create a system image			
Where do you want to sav	the backup?		
A system image is a copy of the dr drives. A system image can be use stops working; however, you can't	es required for Win to restore your con hoose individual ite	dows to run. It can also include additic puter if your hard drive or computer e ms to restore.	onal ever
On a hard disk			
Backup (E:) 465.60 GB free		~	
On one or more DVDs		_	
WD Drive (D:)		~	
On a network location			
		Select	
		Next	Cance

- 5. Verificare le impostazioni di backup e fare clic su Avvia backup.
- 6. Alla richiesta "Creare un disco di ripristino del sistema?", fare clic su No.
- 7. Fare clic su **Chiudi** e rimuovere in modo sicuro l'SSD da Summit. Etichettare l'unità come Ripristino Summit e conservarla in un luogo sicuro.

Ripristino del sistema

In caso di errore del sistema operativo, utilizzare l'unità di ripristino e l'immagine di sistema per ripristinare il sistema.

Materiali richiesti

- Unità di ripristino USB
- SSD con l'immagine di sistema
- Per ripristinare il sistema utilizzando l'unità di ripristino e l'immagine di sistema
- 1. Spegnere lo spettrometro Summit.
- 2. Collegare una tastiera esterna a una delle porte USB presenti sullo spettrometro Summit.
- 3. Collegare l'unità di ripristino USB a una delle porte USB.
- 4. Accendere il sistema.
- 5. Alla comparsa della schermata iniziale Thermo Scientific, toccare "**Esc**" diverse volte sulla tastiera per entrare nel BIOS del sistema.

Se una finestra di dialogo mostra "Quit Without Saving?" (Uscire senza salvare?" utilizzare i tasti di direzione per selezionare No.

- 6. Con i tasti di direzione passare alla scheda Save & Exit (Salva ed esci).
- 7. Passare alla sezione Boot Override (Ignora avvio) e selezionare l'unità flash di ripristino. Premere **Invio** per avviare l'immagine di ripristino.



- 8. Alla comparsa della schermata "Scegli il layout di tastiera", rimuovere l'unità flash di ripristino dalla porta USB e inserire l'SSD dell'immagine di sistema.
- 9. Selezionare il layout di tastiera preferito.
- 10. Nella schermata Scegli un'opzione, selezionare Risoluzione dei problemi.

Iniziare con lo spettrometro Nicolet Summit OA

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

- 11. Selezionare Opzioni avanzate.
- 12. In Opzioni avanzate, selezionare Ripristino immagine di sistema.

	System Restore		Startup Repair
	Use a restore point recorded on your PC to restore Windows	< \$ >	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 file	♦	Startup Settings Change Windows startup behavior

- 13. Selezionare Windows 10 come sistema operativo di destinazione.
- 14. Nella finestra "Selezionare un backup dell'immagine di sistema", utilizzare l'immagine di sistema più recente disponibile oppure scegliere di utilizzare un'altra immagine. Fare clic sul pulsante **Avanti**.

🍋 Re-image your computer			×
	Select a syst	em image backup	
	This computer will b Everything on this information in the s	e restored using the system image. computer will be replaced with the ystem image.	
	Troubleshooting inf http://go.microsofi	ormation for BMR: com/fwlink/p/?LinkId=225039	
		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	
		< <u>B</u> ack <u>N</u> ext > Can	cel

Iniziare con lo spettrometro Nicolet Summit OA

15. Fare clic sul pulsante **Avanti** per mantenere nello stato predefinito ulteriori opzioni di ripristino e fare clic su **Fine** per ripristinare l'immagine di sistema. Fare clic su **Sì** per confermare.

Il ripristino richiede circa 15 minuti e al termine il sistema si riavvia automaticamente.

Riacceso il sistema, rimuovere in modo sicuro l'SSD dallo spettrometro e conservare l'SSD con l'immagine di sistema assieme all'unità di ripristino in un luogo sicuro.

Per informazioni sul ripristino dell'immagine di sistema, vedere la guida in linea di Microsoft.

Backup dei dati di sistema

Oltre a creare supporti di ripristino e un'immagine di sistema, eseguire regolarmente il backup dei dati. Windows 10 offre diverse opzioni per l'esecuzione del backup dei file, incluso lo strumento Cronologia file. Spetta al personale IT dell'organizzazione stabilire quali siano i metodi di backup più adatti alle esigenze dell'organizzazione stessa.

Installazione di un adattatore USB wireless

É possibile collegare lo spettrometro Nicolet Summit OA a una rete wireless utilizzando l'adattatore USB wireless opzionale. Se l'adattatore è stato acquistato con lo spettrometro, è fornito preinstallato.

Per installare l'adattatore USB wireless, inserirlo in una porta USB. Per un'installazione permanente, installare l'adattatore in una porta sotto il coperchio dello strumento.

Per installare un adattatore USB wireless

1. Inserire l'adattatore in una porta USB disponibile.



2. Il sistema riconoscerà l'adattatore automaticamente. Per informazioni su come effettuare il collegamento a una rete utilizzando l'adattatore, vedere il supporto Microsoft.

Per eseguire l'installazione utilizzando una porta USB sotto il coperchio dello strumento

Questa installazione richiede la rimozione della piastra del cristallo e del coperchio dello strumento.

- 1. Seguire le istruzioni contenute in <u>"Riposizionamento del coperchio"</u> per rimuovere la piastra del cristallo e il coperchio dello strumento.
- 2. Con il coperchio rimosso, inserire l'adattatore in una porta USB disponibile.



3. Riposizionare il coperchio dello strumento e la piastra del cristallo.

Lo strumento è pronto per il collegamento a una rete wireless.

Informazioni sulla barra luminosa dello spettrometro

Lo spettrometro Nicolet Summit OA di Thermo Scientific è dotato di una pratica barra luminosa che offre una rapida e immediata indicazione visiva della procedura di test Pass/Fail, per la qualità del prodotto e lo stato dello strumento.

Le tabelle riportate di seguito descrivono tutti i segnali emessi dalla barra luminosa.

RISCALDAMENTO IN CORSO

Segnale	Stato del sistema	Descrizione
Verde ciclico	Sistema acceso e in fase di riscaldamento	La luce verde si sposta ripetutamente da sinistra a destra. Il tempo necessario per il riscaldamento è di circa 2 minuti.

ACQUISIZIONE TRADIZIONALE

Segnale	Stato del sistema	Descrizione
Verde costante	Pronto all'uso	Tutta la barra luminosa mostra una luce verde costante
Cometa verde	Acquisizione dei dati in corso	Una piccola sezione della luce verde rimbalza avanti e indietro

ACQUISIZIONE SMART

Segnale	Stato del sistema	Descrizione
Pulsata blu	Acquisizione dello Smart background	Tutta la barra luminosa emette una luce pulsata blu
Cometa blu	Acquisizione dei dati del campione in corso (iniziata dall'utente)	Una piccola sezione della luce blu rimbalza avanti e indietro

RISULTATI DELL'ANALISI

Segnale	Stato del sistema	Descrizione
% riempimento verde	Valore di corrispondenza o risultato del QCheck sopra la soglia	Riempimento verde per riflettere il valore corrispondente. Ad esempio il riempimento di circa il 90% indica un valore di corrispondenza pari a 90.
% riempimento arancione	Valore di corrispondenza o risultato del QCheck sotto la soglia	Riempimento arancione per riflettere il valore di corrispondenza. Ad esempio il riempimento di circa il 30% indica un valore di corrispondenza pari a 30.

DIAGNOSTICA

Segnale	Stato del sistema	Descrizione
Rossa lampeggiante	Errore di sistema	Tutta la barra luminosa lampeggia in rosso. Per una descrizione dell'errore, vedere Stato del sistema nel software OMNIC Paradigm per una descrizione dell'errore.

Visualizzazione dei dati sulla app OMNIC Anywhere

OMNIC Anywhere è un'applicazione basata sul cloud che consente di visualizzare, esplorare o condividere dati da un PC, un computer Apple o un dispositivo Android o iOS collegato.



Figura 2-1: OMNIC Anywhere nel web browser

Con uno spettrometro Nicolet Summit OA e un account Connect gratuito, potete analizzare i campioni in aula o in laboratorio, caricare i dati sul vostro account Connect e visualizzare, esplorare o condividere i dati su un altro dispositivo del vostro dominio o della vostra area di lavoro.

Per istruzioni su come creare un account Connect e visualizzare i dati su OMNIC, potete visitare "Spettroscopia FTIR abilitata per il cloud". [Pagina lasciata intenzionalmente vuota]

Capitolo 3

Opzioni e accessori

Nicolet Summit OA può essere utilizzato con il kit di spurgo opzionale.

Installazione e manutenzione di un kit di spurgo

Lo spurgo dello spettrometro con aria secca o azoto protegge i componenti interni dall'umidità e altri contaminanti ambientali. L'installazione del kit di spurgo richiede l'assemblaggio di valvole e dispositivi di regolazione, la sostituzione della cartuccia di essiccante, il collegamento del gas di spurgo e l'impostazione della pressione e della portata.

AVVERTENZA



Pericolo di esplosione

Non utilizzare mai un gas infiammabile, combustibile o tossico per spurgare lo strumento. Il gas di spurgo deve essere privo di olio o altri materiali reattivi. Il calore emesso dalla sorgente o dall'assorbimento del laser potrebbe provocare l'accensione di gas infiammabili o di materiali reattivi contenuti nel gas di spurgo. Utilizzare solo aria secca o azoto per spurgare lo strumento.

AVVISO

È consigliabile mantenere sempre la sigillatura e l'essiccante e/o spurgare lo strumento. La garanzia non copre danni derivanti dalla mancata corretta essiccazione o spurgo dello strumento. In caso di domande in merito, rivolgersi alla nostra azienda.

CONFIGURAZIONE RICHIESTA

Prima di installare il kit di spurgo, occorre avere una sorgente di aria secca o azoto compatibile con le specifiche di Thermo Scientific per lo spettrometro Nicolet[™] Summit. Per dettagli, consultare la guida al sito e alla sicurezza dello strumento.

Nota Per ottenere i migliori risultati, seccare il gas di spurgo a un punto di rugiada di -70 °C o inferiore.

MATERIALI RICHIESTI

Oltre al kit di spurgo, occorrono i seguenti strumenti e materiali:

- Chiave a estremità aperta da 3/4"
- Chiave a estremità aperta da 11/16"
- Cacciavite Phillips di misura 2

- Chiave per viti a testa esagonale da 3/32"
- Chiave per viti a testa esagonale da 7/64"
- Nastro sigillante per filettature ("nastro da idraulico" o "Teflon")

Per installare un kit di spurgo

- 1. Collegare il gruppo idraulico alla sorgente del gas di spurgo.
 - a. Installare la valvola regolatrice e un raccordo maschio da 1/4" o femmina da 3/8" sulla sorgente del gas di spurgo (scegliere una valvola e raccordi adeguati per la sorgente del gas di spurgo).



b. In caso di utilizzo di un raccordo maschio da 1/4", procedere alla fase successiva.

In caso di utilizzo di un raccordo femmina da 3/8" sulla sorgente del gas di spurgo, installare il raccordo riduttore da 3/8" a 1/4" contenuto nel kit di spurgo. Prima di installare il raccordo riduttore, avvolgerlo con nastro per filettature e una chiave a estremità aperta da 11/16", stringere la connessione.



c. Avvolgere il raccordo riduttore o il raccordo maschio da 1/4" con il nastro quindi installare l'attacco a pressione. Utilizzare una chiave a estremità aperta da 3/4" per serrare la connessione.



- d. Fare scattare saldamente l'ingresso maschio del gruppo idraulico a parete nell'accessorio a sgancio rapido.
- 2. Spegnere lo spettrometro.
 - a. Per spegnere lo spettrometro, premere il pulsante di alimentazione.
 - b. Scollegare il cavo di alimentazione.
- 3. "Rimozione della piastra del cristallo"
- 4. Rimuovere il coperchio dello strumento.
 - a. Ribaltare con cura lo strumento all'indietro appoggiandolo sulle alette del dissipatore di calore.
 - b. Utilizzando la chiave per viti a testa esagonale da 7/64, allentare le quattro viti che fissano il coperchio alla base dello strumento.



- c. Riportare con cautela lo strumento sui propri piedi. Quando si ribalta nuovamente lo strumento in posizione eretta, le quattro viti cadranno. Fare attenzione a non perderle.
- d. Sollevare il coperchio dritto fuori dalla base.



5. Rimuovere la cartuccia di essiccante.

a. Utilizzare il cacciavite per viti Phillips di misura 2 per allentare le due viti prigioniere dalla cartuccia di essiccante e sollevare verticalmente la cartuccia per estrarla dallo strumento.



b. Rimuovere e scartare l'O-ring grande.



Nota Prestare attenzione per non far cadere l'O-ring nel vano dell'essiccante.

- 6. Aprire il kit di spurgo e la confezione sigillata di essiccante.
- 7. Estrarre l'O-ring dalla confezione e installarlo sulla piastra base dello strumento (vedere immagine precedente). Premere (non ruotare) in posizione il nuovo O-ring in modo da inserirlo correttamente nella scanalatura.

8. Togliere dalla confezione la cartuccia di essiccante con i raccordi di spurgo, inserire la cartuccia nel vano dell'essiccante, assicurarsi del suo corretto inserimento sopra l'O-ring e fissarla mediante le due viti.

Prendere nota dell'orientamento corretto della cartuccia. Quando la cartuccia è installata, deve essere possibile leggere la sua etichetta dalla parte anteriore dello strumento.



- 9. Indirizzare la linea di spurgo attraverso l'apertura dell'ingresso di spurgo sul retro dello strumento.
 - a. Abbassare con cautela il coperchio sopra la base indirizzando al contempo il tubo di spurgo nella fessura del coperchio.



- b. Abbassare completamente il coperchio. Quando il coperchio è abbassato in posizione, il tubo di spurgo è intrappolato nella fessura fra il coperchio e il dissipatore di calore.
- c. Con il tubo di spurgo spinto di lato, ribaltare lo strumento all'indietro, appoggiandolo sulle alette del dissipatore di calore.
- d. Inserire le quattro viti precedentemente rimosse. Serrare tutte le viti.
- e. Riportare nuovamente lo strumento sui propri piedi.
- f. Spingere lo spezzone di tubo in eccesso nello strumento o estrarlo secondo necessità.



10. "Riposizionamento della piastra del cristallo".

- 11. Collegare il giunto di flusso dal gruppo idraulico a parete (indicato con "allo spettrometro") al tubo di spurgo.
- 12. Impostare i comandi del gas di spurgo.
 - a. Aprire la valvola regolatrice principale e impostare i comandi del gas di spurgo come sotto indicato (vedere Impostazione dei comandi del gas di spurgo per dettagli).

Hardware	Pressione (psig)	Portata (scfh)
Spettrometro Summit	5	1

- b. Collegare il cavo di alimentazione allo strumento e accendere l'alimentazione.
- c. Attendere per 30-60 minuti lo spurgo completo dello strumento.

Impostazione dei comandi del gas di spurgo

Impostando correttamente i comandi del gas di spurgo è possibile proteggere lo spettrometro dall'umidità senza introdurre vibrazioni.

Per ottenere i migliori risultati, seccare il gas di spurgo a una temperatura di condensazione di - 70 °C o inferiore.



- 1. Aprire la valvola principale per avviare il flusso del gas di spurgo attraverso il regolatore.
- 2. Regolare il regolatore di pressione per lo spettrometro finché il misuratore non indica una pressione di 5 psig (34 kPa).
- 3. Impostare il flussimetro per lo spettrometro su 1 scfh (0,47 l/min).

AVVISO

Portate superiori ai valori raccomandati possono causare una vibrazione che può influire sulla qualità dei dati.

Controllo e sostituzione del filtro del gas di spurgo

Sostituire il filtro di spurgo quando è giallo, o altrimenti scolorito, o se è contaminato da detriti o particelle estranee.

4.

AVVISO

È consigliabile mantenere sempre sigillato ed essiccato e/o spurgare lo strumento. I danni all'apparecchiatura dovuti al mancato mantenimento della sigillatura e dell'essiccazione e/o spurgo non sono inclusi nella garanzia. In caso di domande in merito, rivolgersi alla nostra azienda.

Figura 3-1: Il filtro di spurgo si trova all'interno del bicchierino in plastica sotto il manometro



Nota Per ordinare componenti, contattare il Supporto tecnico.

Per sostituire il filtro del gas di spurgo

- 1. Chiudere il gas di spurgo dalla valvola principale. Non capovolgere il flussimetro o il regolatore di pressione.
- 2. Rimuovere il bicchierino in plastica che alloggia il filtro e quindi rimuovere il filtro (svitandoli entrambi manualmente).



- 3. Installare il nuovo filtro e reinstallare il bicchierino.
- 4. Aprire la valvola principale per attivare il flusso di spurgo allo strumento e verificare che la portata del gas di spurgo sia appropriata.

[Pagina lasciata intenzionalmente vuota]

Capitolo 4

Manutenzione

Lo spettrometro Nicolet Summit OA FTIR richiede un modesto mantenimento e nessuna manutenzione quotidiana. Tuttavia la sua cura di base garantisce costantemente prestazioni di massimo livello.

Pianificazione della manutenzione per lo spettrometro Summit	
Allineamento delle ottiche dello spettrometro	
Calibrazione della frequenza del laser	
Pulizia dello strumento e del cristallo	
Rimozione e riposizionamento della piastra del cristallo	
Riposizionamento del coperchio	54
Sostituzione dell'essiccante	
Sostituzione della sorgente	

Pianificazione della manutenzione per lo spettrometro Summit

Lo spettrometro Nicolet Summit OA FTIR di Thermo Scientific richiede un modesto mantenimento e nessuna manutenzione quotidiana. Tuttavia la sua cura di base garantisce costantemente prestazioni di massimo livello. Seguire le linee guida fornite in questa sezione. Per ottenere prestazioni ottimali, lasciare acceso lo spettrometro.

AVVISO

L'elettricità statica può danneggiare in modo permanente componenti critici dello spettrometro. Per contribuire a evitare un danno di questo tipo, seguire queste raccomandazioni:

- Prima di staccare l'alimentazione, scaricare l'elettricità statica che potrebbe essersi accumulata, toccando la base metallica dello spettrometro.
- Lasciare i componenti di ricambio nei rispettivi imballaggi di protezione fino al momento di installarli nello strumento.

Manutenzione settimanale

VERIFICA DELLE PRESTAZIONI DELLO SPETTROMETRO

Il software OMNIC Paradigm include un workflow per la verifica delle prestazioni per lo spettrometro. Il workflow per la verifica delle prestazioni (PV) esegue una serie di test standard volti a verificare il funzionamento dello strumento e garantire l'accuratezza dei dati.

È consigliabile eseguire il workflow PV almeno una volta alla settimana.

PULIZIA DELLO SPETTROMETRO

Lo spettrometro può essere pulito solo in conformità alle raccomandazioni. Vedere <u>"Pulizia dello</u> spettrometro".

Manutenzione mensile

CONTROLLO DELL'INDICATORE DI UMIDITÀ

Le ottiche racchiuse nello spettrometro possono venire facilmente danneggiate da un'eccessiva umidità dell'aria. Lo spettrometro è sigillato e i suoi componenti protetti da due contenitori di essiccante in grado di assorbire l'umidità. Monitorare l'umidità dello strumento nel software OMNIC Paradigm.

AVVISO

È consigliabile mantenere sempre lo spettrometro sigillato ed essiccato oppure sigillato e spurgato con aria secca o azoto. I danni all'apparecchiatura dovuti al mancato mantenimento della sigillatura e dell'essiccazione e/o spurgo non sono inclusi nella garanzia. In caso di domande in merito, rivolgersi alla nostra azienda.

CONTROLLO DEL FILTRO DEL GAS DI SPURGO

Se lo spettrometro è spurgato con azoto o aria secca, controllare il filtro di spurgo almeno una volta al mese.

Sostituire il filtro quando è giallo, o altrimenti scolorito, o se è contaminato da detriti o particelle estranee. Per ulteriori informazioni, vedere <u>"Controllo e sostituzione del filtro del gas di spurgo"</u>.

Figura 4-1: Il filtro di spurgo si trova all'interno del bicchierino in plastica sotto il manometro



Allineamento delle ottiche dello spettrometro

Allineare le ottiche dello spettrometro qualora il workflow di verifica delle prestazioni Nicolet Summit OA ne indichi il mancato superamento. L'allineamento dello spettrometro permette di ottimizzare l'energia che raggiunge il rivelatore e di massimizzarne il segnale. Se il test delle prestazioni non viene superato, è necessario allineare le ottiche dello spettrometro, calibrare la frequenza del laser e rieseguire il test. Se il test non viene superato nemmeno in questo caso, contattare il rappresentante del supporto tecnico locale per chiedere assistenza.

- Per allineare le ottiche dello spettrometro (interfaccia touchscreen)
- 1. Nella schermata iniziale, selezionare l'icona Diagnostica [200] per aprire la vista omonima.
- 2. Aprire la scheda Allinea e selezionare Allinea.

Il sistema allinea automaticamente lo spettrometro e indica quando l'operazione è stata completata.

- Per allineare le ottiche dello spettrometro (interfaccia operatore)
- 1. Spostarsi nell'area Amministratore, selezionare l'icona Diagnostica [
- 2. Aprire la scheda Allinea e selezionare Allinea.

Il sistema allinea automaticamente lo spettrometro e indica quando l'operazione è stata completata.

- Per allineare le ottiche dello spettrometro (interfaccia desktop)
- 1. Selezionare Acquisisci dati > Diagnostica > Allinea spettrometro.
- 2. Nella finestra di dialogo Allinea spettrometro, fare clic su Avvio.

Il sistema allinea automaticamente lo spettrometro e indica quando l'operazione è stata completata.

Calibrazione della frequenza del laser

Calibrare la frequenza del laser se il workflow di verifica delle prestazioni Nicolet Summit OA indica il mancato superamento del test. Se un test non viene superato, è necessario prima allineare lo spettrometro, quindi calibrare la frequenza del laser e rieseguire il test. Se il test non viene superato nemmeno in questo caso, contattare il rappresentante del supporto tecnico locale per chiedere assistenza.

Per calibrare il laser (interfaccia touchscreen)

- 1. Nella schermata iniziale, selezionare l'icona Diagnostica [
- 2. Aprire la scheda Calibrazione laser e selezionare Calibra per cominciare.

Completata la calibrazione del laser, un messaggio indica se la frequenza del laser è stata modificata e visualizza la nuova frequenza.

Per calibrare il laser (interfaccia operatore)

- 1. Spostarsi nell'area Amministratore, selezionare l'icona Diagnostica [
- 2. Aprire la scheda Calibrazione laser e selezionare Calibra per cominciare.

Completata la calibrazione del laser, un messaggio indica se la frequenza del laser è stata modificata e visualizza la nuova frequenza.

Per calibrare il laser (interfaccia desktop)

- 1. Selezionare Acquisisci dati > Diagnostica > Calibrazione laser.
- 2. Nella finestra di dialogo Calibrazione laser, fare clic su Avvio per calibrare il laser.

Quando il sistema ha terminato la calibrazione del laser, un messaggio indica se la frequenza è stata modificata e visualizza la nuova frequenza.

Pulizia dello strumento e del cristallo

Prima di pulire lo spettrometro Nicolet Summit OA di Thermo Scientific, spegnerlo e staccare l'alimentazione.

Pulizia dello spettrometro

Pulire delicatamente la superficie esterna dello strumento utilizzando un panno morbido pulito e un solvente delicato come Windex. Spruzzare il solvente sul panno e pulire solo la superficie esterna dipinta; evitare il contatto con i componenti elettronici o la parte posteriore dello strumento.

Sull'involucro dei componenti elettronici sul retro dello strumento si può accumulare polvere che, interferendo potenzialmente con la dissipazione del calore, può ridurre la vita utile dei componenti elettronici.

Rimuovere la polvere dalla parte posteriore dello strumento utilizzando un soffio di aria compressa. Non utilizzare liquidi per rimuovere la polvere dalla parte posteriore dello spettrometro.

ATTENZIONE



Pericolo di scosse elettriche.Per evitare il pericolo di scosse elettriche, occorre impedire la penetrazione di liquidi nell'alimentatore o sul retro dello strumento.

AVVISO

Non utilizzare detergenti aggressivi, solventi, sostanze chimiche o abrasive che potrebbero danneggiare la finitura.

Pulizia della piastra del cristallo ATR

Utilizzare un panno o una salvietta non abrasivi puliti e asciutti per pulire il cristallo ATR.

AVVERTENZA



Pericolo di inalazione e ingestione di sostanze tossiche.

Il seleniuro di zinco (ZnSe) è tossico. Fare riferimento alla Scheda di dati di sicurezza ZnSe del produttore all'indirizzo www.specac.com per la manipolazione e i controlli dell'esposizione.

Indossare guanti di protezione per pulire o sostituire la piastra del cristallo.

- 1. Utilizzare un panno o una salvietta puliti per assorbire il liquido dal cristallo. Gettare la salvietta usata.
- 2. Con una pipetta di trasferimento, applicare alcune gocce di eptano sul cristallo e sulla zona circostante.
- 3. Con una salvietta nuova pulita, strofinare l'intera lunghezza del cristallo per rimuovere interamente i liquidi. Strofinare la zona circostante la vaschetta del cristallo.



4. Ripetere secondo necessità, terminando sempre con un panno o una salvietta puliti e asciutti.

Rimozione e riposizionamento della piastra del cristallo

La piastra del cristallo deve essere sostituita ogni 12-18 mesi o secondo necessità.

AVVERTENZA



Pericolo di inalazione e ingestione di sostanze tossiche.

Il seleniuro di zinco (ZnSe) è tossico. Fare riferimento alla Scheda di dati di sicurezza ZnSe del produttore all'indirizzo www.specac.com per la manipolazione e i controlli dell'esposizione.

Indossare guanti di protezione per pulire o sostituire la piastra del cristallo.

AVVISO

Fare attenzione per evitare che cadano oggetti all'interno dello strumento quando il coperchio è stato rimosso.

MATERIALI RICHIESTI

• Chiave per viti a testa esagonale da 3/32"

Rimozione della piastra del cristallo

- 1. Rimuovere il coperchio antipolvere dal cristallo.
- 2. Utilizzando la chiave per viti a testa esagonale da 3/32" allentare completamente le tre viti che fissano la piastra del cristallo allo strumento.



- 3. Sollevare la piastra del cristallo dritta fuori dallo strumento.
- 4. Utilizzare un panno pulito e asciutto per pulire la superficie circostante la flangia, sotto la zona di appoggio della piastra.

Riposizionamento della piastra del cristallo

- 1. Posizionare la nuova piastra sopra la flangia.
- 2. Applicare le rondelle sulle viti e inserire le viti nei tre fori. Le viti devono essere inserite dritte nei fori.
- 3. Premendo sulla piastra del cristallo, stringere le due viti ai lati della piastra, quindi serrare la terza vite più vicina alla parte posteriore dello strumento. Stringere tutte le viti serrandole a mano.

Se la vite ruota liberamente, estrarla e reinserirla per correggere l'allineamento.

Riposizionamento del coperchio

STRUMENTI RICHIESTI

Prima di cominciare, occorre avere a disposizione i seguenti strumenti:

- Chiave per viti a testa esagonale da 7/64"
- Chiave per viti a testa esagonale da 3/32"
- Per riposizionare il coperchio dello spettrometro
- 1. Spegnere lo spettrometro.
 - a. Per spegnere lo spettrometro, premere il pulsante di alimentazione.
 - b. Scollegare il cavo di alimentazione.
 - c. Disattivare lo spurgo, ove opportuno, e staccare la linea di spurgo collegata allo strumento.
- 2. Rimuovere la piastra del cristallo.
 - a. Rimuovere il coperchio antipolvere dal cristallo.
 - b. Allentare completamente le tre viti che fissano la piastra del cristallo allo strumento.



- c. Sollevare la piastra del cristallo dritta fuori dallo strumento.
- d. Utilizzare un panno pulito e asciutto per pulire la superficie circostante la flangia, sotto la zona di appoggio della piastra.
- 3. Rimuovere il coperchio dello strumento.
 - a. Ribaltare con cura lo strumento all'indietro appoggiandolo sulle alette del dissipatore di calore.
 - b. Utilizzando la chiave per viti a testa esagonale da 7/64, allentare le quattro viti che fissano il coperchio alla base dello strumento.



c. Riportare con cautela lo strumento sui propri piedi. Quando si ribalta nuovamente lo strumento in posizione eretta, le quattro viti cadranno. Fare attenzione a non perderle.

d. Sollevare il coperchio dritto fuori dalla base.



- 4. Riposizionare il coperchio dello strumento.
 - a. Portare con cautela il coperchio in posizione sulla base dello strumento.
 - b. Ribaltare lo strumento all'indietro appoggiandolo sulle alette del dissipatore di calore.
 - c. Inserire le quattro viti precedentemente rimosse. Serrare tutte le viti.
 - d. Riportare nuovamente lo strumento sui propri piedi.
- 5. Riposizionare la piastra del cristallo.
 - a. Posizionare la piastra sopra la flangia.
 - b. Applicare le rondelle sulle viti e inserire le viti nei tre fori. Le viti devono essere inserite dritte nei fori.
 - c. Premendo sulla piastra del cristallo, stringere le due viti ai lati della piastra, quindi serrare la terza vite più vicina alla parte posteriore dello strumento. Stringere tutte le viti serrandole a mano.

Se la vite ruota liberamente, estrarla e reinserirla per correggere l'allineamento.

6. Accendere lo spettrometro.

- a. Collegare il cavo di alimentazione allo strumento e accendere l'alimentazione.
- b. Ricollegare la linea di spurgo allo strumento, se pertinente, e attivare lo spurgo.
- 7. Verificare le prestazioni dello spettrometro.
 - a. Allineamento delle ottiche dello spettrometro.
 - Dal dashboard del software OMNIC Paradigm andare a Acquisisci dati > Diagnostica > Allinea spettrometro.
 - ii. Selezionare **Avvio** per iniziare l'allineamento. La schermata visualizza "Allineamento completato" quando l'allineamento è terminato con successo.
 - b. Calibrare il laser.
 - i. Andare a Acquisisci dati > Diagnostica > Calibrazione laser.
 - ii. Selezionare Avvio per iniziare la calibrazione.

Il completamento della calibrazione laser può richiedere diversi minuti. Terminata la calibrazione, il display visualizza "Misura completata" e indica le modifiche della frequenza laser.

Lo strumento impiega 12 ore per raggiungere l'equilibrio termico. Per conseguire i migliori risultati, rieseguire l'allineamento e la calibrazione dopo le prime 12 ore di funzionamento.

- c. Eseguire il **workflow di verifica delle prestazioni Nicolet Summit OA**. Il workflow di verifica delle prestazioni utilizza materiale di riferimento per eseguire una serie di prove standard per verificare il funzionamento dello strumento e garantire l'accuratezza dei dati.
 - i. Ritornare al Dashboard.
 - ii. Scorrere verso il riquadro Workflow e selezionare Verifica delle prestazioni Nicolet Summit OA.
 - iii. Fare clic con il tasto destro sul workflow e selezionare **Esegui** per avviare il workflow.

Sostituzione dell'essiccante

Monitorare l'umidità dello strumento utilizzando il software OMNIC Paradigm. Sostituire l'essiccante secondo necessità o attenendosi ai criteri stabiliti dalla propria organizzazione.

Per sostituire l'essiccante, rimuovere il coperchio dello strumento, la cartuccia dell'essiccante e sostituire i contenitori di essiccante e l'O-ring.

STRUMENTI RICHIESTI

Prima di cominciare, occorre avere a disposizione i seguenti strumenti:

- Chiave per viti a testa esagonale da 7/64"
- Kit di sostituzione dell'essiccante
- Chiave per viti a testa esagonale da 3/32"
- Cacciavite Phillips di misura 2
- Guanti, protezioni per le dita o salvietta da laboratorio (per manipolare l'indicatore di umidità)

AVVERTENZA



Pericolo di esplosione.

Prima di rimuovere il coperchio dello strumento e aprire il vano dell'essiccante, spegnere lo strumento, scollegare il cavo di alimentazione e rimuovere i campioni dal sistema.

L'ingresso di liquidi o gas infiammabili nel vano dell'essiccante rappresenta un pericolo di esplosione. In tale evenienza, contattare immediatamente la nostra azienda e non alimentare elettricamente lo strumento finché la condizione non sia stata risolta.

AVVISO

Quando il coperchio dell'essiccante viene rimosso, verificare che non cada nulla all'interno dello strumento.

AVVERTENZA



Sostituire l'essiccante solo con ricambi forniti dalla nostra azienda.

Thermo Scientific

Per sostituire l'essiccante

- 1. Spegnere lo spettrometro.
 - a. Per spegnere lo spettrometro, premere il pulsante di alimentazione.
 - b. Scollegare il cavo di alimentazione.
 - c. Disattivare lo spurgo, ove opportuno, e staccare la linea di spurgo collegata allo strumento.

Nota È possibile utilizzare una cartuccia di essiccante (con o senza il connettore di spurgo) per preservare l'umidità richiesta. Le immagini di seguito riportate mostrano la versione della cartuccia di essiccante non spurgata. Le istruzioni per la sostituzione dei contenitori di essiccante sono le stesse per entrambe le cartucce.

- 2. "Rimozione della piastra del cristallo"
- 3. Rimuovere il coperchio dello strumento.
 - a. Ribaltare con cura lo strumento all'indietro appoggiandolo sulle alette del dissipatore di calore.
 - b. Utilizzando la chiave per viti a testa esagonale da 7/64, allentare le quattro viti che fissano il coperchio alla base dello strumento.



c. Riportare con cautela lo strumento sui propri piedi. Quando si ribalta nuovamente lo strumento in posizione eretta, le quattro viti cadranno. Fare attenzione a non perderle.

d. Sollevare il coperchio dritto fuori dalla base.



- 4. Rimuovere la cartuccia di essiccante.
 - a. Utilizzare il cacciavite per viti Phillips di misura 2 per allentare le due viti prigioniere dalla cartuccia di essiccante e sollevare verticalmente la cartuccia per estrarla dallo strumento.



b. Rimuovere e scartare l'O-ring grande.



Nota Prestare attenzione per non far cadere l'O-ring nel vano dell'essiccante.

5. Rimuovere le cartucce dell'essiccante e sostituire l'indicatore di umidità.



Il contenuto dei contenitori di essiccante (setaccio molecolare alluminosilicato metallico idratato) può essere nocivo in caso di ingestione. Accertarsi che i contenitori di essiccante saturo vengano correttamente smaltiti.

a. Utilizzare la chiave esagonale da 0,05" per allentare (circa due giri) le otto viti che tengono i due contenitori di essiccante saturo sulla relativa cartuccia. Rimuovere i contenitori.



b. Capovolgere la cartuccia di essiccante e staccare il vecchio indicatore di umidità dalla finestra. Gettare l'indicatore usato.

AVVISO

Indossare sempre guanti da laboratorio o protezioni per le dita, o utilizzare una salvietta da laboratorio, per maneggiare un indicatore di umidità nuovo. Il grasso o l'umidità della cute possono scolorire l'indicatore.

- 6. Installare i nuovi contenitori di essiccante.
 - a. Aprire il pacchetto sigillato del nuovo essiccante (deve essere sigillato per preservare l'essiccante) e rimuovere il nuovo indicatore di umidità.
 - b. Capovolgere la cartuccia e premere l'indicatore blu sulla finestra in modo che i bordi piatti dell'indicatore siano allineati ai bordi della cartuccia. Quando la cartuccia viene raddrizzata il testo deve essere visibile attraverso la finestra.

AVVISO

Accertarsi che l'indicatore sia premuto saldamente contro la finestra in modo che non si allenti nel vano dell'essiccante durante la reinstallazione della cartuccia.

- c. Inserire i contenitori nel canale della cartuccia di essiccante e stringere le otto viti (circa 2 giri) sino a portarli al livello del canale della cartuccia.
- 7. Inserire la cartuccia di essiccante.
 - a. Estrarre il nuovo O-ring dal pacchetto e installarlo sulla piastra base dello strumento.

Premere, anziché ruotare in posizione, il nuovo O-ring in modo da inserirlo correttamente nella scanalatura.

b. Inserire la cartuccia di essiccante nello spettrometro in modo che il testo riportato sull'indicatore di umidità sia leggibile quando ci si trova davanti allo strumento.

Accertarsi che la cartuccia sia correttamente inserita in sede sopra l'O-ring quindi utilizzare il cacciavite per viti Phillips di misura 2 per fissare le due viti prigioniere.

- 8. Riposizionare il coperchio dello strumento.
 - a. Portare con cautela il coperchio in posizione sulla base dello strumento.
 - b. Ribaltare lo strumento all'indietro appoggiandolo sulle alette del dissipatore di calore.
 - c. Inserire le quattro viti precedentemente rimosse. Serrare tutte le viti.

- d. Riportare nuovamente lo strumento sui propri piedi.
- e. "Riposizionamento della piastra del cristallo".
- 9. Accendere lo spettrometro.
 - a. Collegare il cavo di alimentazione allo strumento e accendere l'alimentazione.
 - b. Ricollegare la linea di spurgo allo strumento, se pertinente, e attivare lo spurgo.
- 10. Verificare le prestazioni dello spettrometro.
 - a. Allineamento delle ottiche dello spettrometro.
 - i. Dal dashboard del software OMNIC Paradigm andare a Acquisisci dati > Diagnostica > Allinea spettrometro.
 - ii. Selezionare **Avvio** per iniziare l'allineamento. La schermata visualizza "Allineamento completato" quando l'allineamento è terminato con successo.
 - b. Calibrare il laser.
 - i. Andare a Acquisisci dati > Diagnostica > Calibrazione laser.
 - ii. Selezionare Avvio per iniziare la calibrazione.

Il completamento della calibrazione laser può richiedere diversi minuti. Terminata la calibrazione, il display visualizza "Misura completata" e indica le modifiche della frequenza laser.

Lo strumento impiega 12 ore per raggiungere l'equilibrio termico. Per conseguire i migliori risultati, rieseguire l'allineamento e la calibrazione dopo le prime 12 ore di funzionamento.

- c. Eseguire il **workflow di verifica delle prestazioni Nicolet Summit OA**. Il workflow di verifica delle prestazioni utilizza materiale di riferimento per eseguire una serie di prove standard per verificare il funzionamento dello strumento e garantire l'accuratezza dei dati.
 - i. Ritornare al Dashboard.
 - ii. Scorrere verso il riquadro Workflow e selezionare Verifica delle prestazioni Nicolet Summit OA.
 - iii. Fare clic con il tasto destro sul workflow e selezionare **Esegui** per avviare il workflow.

Sostituzione della sorgente

La sorgente IR è facilmente accessibile dal fondo dello spettrometro e per sostituirla non è nemmeno necessario rimuovere il coperchio.

STRUMENTI E MATERIALI RICHIESTI

- Un cacciavite Phillips di misura 1
- Kit di sostituzione della sorgente

ATTENZIONE



Prima di sostituire la sorgente, spegnere lo strumento. Sostituire la sorgente solo con ricambi forniti dalla nostra azienda.

Per sostituire la sorgente

- 1. Spegnere e scollegare lo spettrometro.
 - a. Spegnere lo strumento e staccare il cavo di alimentazione.
 - b. Scollegare dallo strumento tutti i cavi, come il cavo Ethernet o i dispositivi USB.
 - c. Se lo spettrometro è spurgato, staccare le linee di spurgo. Per ulteriori informazioni, vedere "Installazione e manutenzione di un kit di spurgo".

ATTENZIONE



d.

Pericolo di ustioni.

La sorgente diventa molto calda durante l'uso dello strumento. Attendere almeno 10 minuti dopo lo spegnimento dello strumento prima di eseguire la fase successiva.

- 2. Rimuovere la sorgente.
 - a. Con un cacciavite Phillips misura 1 allentare completamente le tre viti prigioniere che tengono in posizione la sorgente.



b. Far fuoriuscire l'involucro. Tenere fermo lo strumento con una mano tirando saldamente l'involucro per estrarre la sorgente dallo strumento.



AVVISO

Non toccare mai un elemento sorgente a mani nude. I grassi cutanei o altri depositi sull'elemento ne riducono la durata. Utilizzare sempre protezioni per mantenere le dita pulite, guanti o una salvietta da laboratorio pulita per maneggiare l'elemento sorgente.

- 3. Inserire la nuova sorgente.
 - a. Accertarsi che le tre viti prigioniere siano diritte quindi inserire lentamente la sorgente nella cavità montandola a filo con lo strumento.
 - b. Stringere le viti per fissare la sorgente e fare scattare l'involucro in posizione di stoccaggio.
- 4. Ricollegare e accendere lo spettrometro.
 - a. Riportare con cautela lo strumento in posizione verticale e ricollegare tutti i cavi rimossi.
 - b. Sostituire l'essiccante. Per istruzioni vedere "Sostituzione dell'essiccante".

Poiché rimuovendo la sorgente le ottiche interne vengono esposte all'atmosfera, l'essiccante deve sempre essere sostituito dopo la sostituzione della sorgente.

AVVERTENZA



Pericolo di scosse elettriche.

Per assicurare una buona messa a terra ed evitare il pericolo di scosse, non usare una presa collegata a una canalizzazione di messa a terra. La terra deve essere un conduttore che non porta corrente, con collegamento a terra nel quadro di distribuzione principale.

- c. Collegare il cavo di alimentazione allo strumento e premere l'interruttore di alimentazione per accendere lo spettrometro.
- 5. Verificare le prestazioni dello spettrometro.
 - a. Allineamento delle ottiche dello spettrometro.
 - i. Dal dashboard del software OMNIC Paradigm andare a Acquisisci dati > Diagnostica > Allinea spettrometro.
 - ii. Selezionare **Avvio** per iniziare l'allineamento. La schermata visualizza "Allineamento completato" quando l'allineamento è terminato con successo.
 - b. Calibrare il laser.
 - i. Andare a Acquisisci dati > Diagnostica > Calibrazione laser.
 - ii. Selezionare Avvio per iniziare la calibrazione.

Il completamento della calibrazione laser può richiedere diversi minuti. Terminata la calibrazione, il display visualizza "Misura completata" e indica le modifiche della frequenza laser.

Lo strumento impiega 12 ore per raggiungere l'equilibrio termico. Per conseguire i migliori risultati, rieseguire l'allineamento e la calibrazione dopo le prime 12 ore di funzionamento.

- c. Eseguire il **workflow di verifica delle prestazioni Nicolet Summit OA**. Il workflow di verifica delle prestazioni utilizza materiale di riferimento per eseguire una serie di prove standard per verificare il funzionamento dello strumento e garantire l'accuratezza dei dati.
 - i. Ritornare al Dashboard.
 - ii. Scorrere verso il riquadro Workflow e selezionare Verifica delle prestazioni Nicolet Sum-

mit OA.

iii. Fare clic con il tasto destro sul workflow e selezionare **Esegui** per avviare il workflow.

[Pagina lasciata intenzionalmente vuota]