Nicolet Summit OA FTIR 분광기





269-346500개정 B2021년 4월

© 2020 Thermo Fisher Scientific Inc. All rights reserved.

기술 지원이 필요한 경우 www.thermofisher.com으로 문의하십시오.

Thermo Fisher Scientific Inc.는 제품 구매 고객에게 제품 작동 시 사용하도록 이 문서를 제 공합니다. 이 문서는 저작권으로 보호되어 있으며, Thermo Fisher Scientific Inc.의 서면 승 인이 있는 경우를 제외하고 이 문서의 전체 또는 일부를 복제하는 행위는 엄격히 금지됩 니다.

이 문서의 내용은 통지 없이 변경될 수 있습니다.이 문서의 모든 기술 정보는 참조용으 로만 제공됩니다.이 문서의 시스템 구성과 사양은 구매자가 받은 이전의 정보를 모두 대 체합니다.

Thermo Fisher Scientific Inc.는 이 문서가 완벽하거나 정확하거나 오류가 없다는 표현 을 하지 않으며, 이 문서의 정보를 적절하게 따랐다 하더라도 이 문서의 사용으로 초래 될 수 있는 오류, 누락, 손상 또는 손실에 대해 일체 그 책임을 지지 않습니다.

이 문서는 Thermo Fisher Scientific Inc.와 구매자 간 판매 계약의 일부가 아닙니다. 이 문서 는 어떤 형태로든 판매 조건을 통제하거나 수정하지 않으며, 판매 조건은 두 문서 간에 상충하는 모든 정보를 통제합니다.

연구용으로만 사용하십시오 이 장비 또는 부속품은 의료 기기가 아니며, 질병의 예방, 진단 또는 치료에 사용하는 용도로 제작되지 않았습니다.



목차

시작	1
사용된 표기법	2
부품 주문	3
연락처	3

Nicolet Summit OA 분광기 시작하기	5
Summit 분광기가 도착하기 전에	6
Summit 분광기와 개인적 안전	
분광기 포장 풀기 및 시작하기	
데이터 백업 및 복구	
USB 무선 어댑터 설치	24
분광계 LightBar 이해	
OMNIC Anywhere 앱으로 데이터 보기	

옵션 및 액세서리	
퍼지 킷 설치 및 유지보수	32

유지보수	41
Summit 분광기의 유지보수 일정	42
분광기 광학장치 정렬	
레이저 주파수 보정	
장비와 크리스탈 청소	46
크리스탈 플레이트 탈거 및 교환	48
덮개 교환	
건조제 교환	54
소스 교환	60

시작

Thermo Scientific Nicolet Summit OA FTIR(Fourier Transform Infrared) 분광기를 사용하면 중간 IR 스펙트럼 범위에서 데이터를 수집하여, 샘플 물질의 화학 분석을 수행할 수 있습니다. 이 시스 템에는 확인 기능, 강력한 소프트웨어 세트 그리고, 데이터 수집을 간편하게 해 주는 다른 많은 기능이 통합되어 있습니다. 옵션 하드웨어를 설치하고, 여러 서비스와 유지보수 절차를 직접 수행할 수 있습니다. 이 문서 또는 제공된 다른 설명서에 사용자에게 필요한 정보가 들어 있습 니다.

장비를 사용하기 전에 시스템과 함께 제공된 안전 설명서를 참조하십시오.

참고 항상 씰과 건조를 유지하거나 장비를 퍼지하는 것이 좋습니다. 씰과 건조를 유지하거 나 퍼징하지 못한 결과로 인한 장비 손상에는 보증이 적용되지 않습니다. 이 요건에 관하여 궁금한 점이 있는 경우 당사에 문의하십시오.

사용된 표기법

안전 주의사항 및 기타 중요한 정보는 다음 형식을 사용합니다.



부품 주문

부품을 주문하려면 당사에 문의하십시오.

장비 또는 부속품을 당시에 수리하기 위해 보내야 하는 경우 배송 요건이나 다른 지침이 필요하 다면 먼저 당사에 전화하거나 이메일을 보내 주세요.

연 락 처

기술 지원이 필요한 경우 www.thermofisher.com으로 문의하십시오.

[별도 공란]

Nicolet Summit OA 분광기 시작 하기

새 Nicolet Summit OA 분광기를 시작하려면 장비가 도착하기 전에 작업장 및 안전 요건을 검토 하고 장비 포장을 풀고 설치한 후 추가적인 부속품을 연결하십시오.

Summit 분광기가 도착하기 전에	6
Summit 분광기와 개인적 안전	
분광기 포장 풀기 및 시작하기	14
데이터 백업 및 복구	
USB 무선 어댑터 설치	24
분광계 LightBar 이해	26
OMNIC Anywhere 앱으로 데이터 보기	

Summit 분광기가 도착하기 전에

데이터의 정확도와 장기적인 성능이 극대화되도록 하려면 도착 전에 적당한 분광기 작업 공간 을 준비하십시오.

이 항목에서는 분광기가 도착하면 할 일을 설명하며, 분광기의 성능에 영향을 미칠 수 있는 환 경/전기적 요소에 대해 소개합니다. 이러한 요건에 대한 더 포괄적인 설명은 Summit 분광기의 작업장 및 안전 정보를 참조하십시오.

분광기 포장 풀기

상자를 개봉하기 전에,장비 도착 즉시 수행해야 할 2가지 중요한 단계가 있습니다.

• 제품 상자의 외부에 손상이 있는지 여부를 확인하십시오.

손상이 있는 경우 당사에 또는 해당하는 현지 판매점에 지침을 문의하십시오.

• 분광기가 실온에 도달하도록 두십시오.

제품 상자 내부의 분광기는 플라스틱 백으로 밀봉되어 건조 상태를 유지하고 있습니다. 백 을 개봉하기 전에 분광기가 실온에 도달하도록 24시간 동안 그대로 두어야 합니다.</u>분광기 가 예열되기 전에 백을 개봉하면 내부 광학 구성부품을 손상시킬 수 있는 응축이 형성되고 영구적인 손상이 발생할 수 있습니다.

장비가 실온에 도달하기 전에 밀봉된 플라스틱 백을 개봉하거나 이동 시 잘못 취급하여 발생한 손상에는 보증이 적용되지 않습니다.

작업 공간 준비

장비가 도착하기 전에, 작업 공간이 분광기를 올바르게 지원할 수 있는지 확인하십시오. 분광 기를 사용하기 위한 충분한 공간을 확보하는 것 외에, 여러 가지 환경 그리고 전기적 요건을 고 려하십시오.

분광기 사이즈

Summit 분광기에 필요한 공간은 비교적 작지만, 열이 환기구에서 분산되고 사용자가 장비의 포 트, 전원, 케이블에 쉽게 접근할 수 있도록 장비 주위에도 여유 공간을 두십시오.

- Summit 장비 무게: 12.6 kg(27.8 파운드)
- 사이즈(너비 x 높이 x 깊이) 34 cm x 24 cm x 32 cm(13.3 인치 x 9.6 인치 x 12.7 인치)

환경 요소

Summit 분광기는 다양한 환경에 사용하도록 설계된 강력한 장비입니다.그러나 최고의 성능을 유지하려면 상대적으로 먼지가 덜하고 습도가 낮은 환경에 보관하십시오.이 분광기는 15°C~ 35°C의 온도에서 안정적으로 작동하지만, 최적의 성능을 위해서는 20°C~22°C로 온도를 유지 해야 합니다.

습기는 장비 내부에 응축을 유발하여, 내부 구성부품이 손상될 수 있습니다. 보통은 다음과 같 은 몇 가지 예방 조치로 장비를 습기로부터 보호할 수 있습니다.

- 장비를 보관 중일 때를 포함하여 장비의 건조제를 유지합니다.
- 급격한 온도 변화를 피합니다.
 - 장비를 냉난방 환기구 또는 큰 창문 근처와 같이 차거나 뜨거운 공기 공급원에 가까이 두지 마십시오.

분광기를 특히 습한 환경에 보관하게 될 경우 퍼지 가스 킷 설치를 고려하십시오.

전기 요건

분광기에 공급되는 전원은 전용 무정전 전원이고 다음 문제가 없어야 합니다.

- 전압강하
- 일시적 스파이크
- 주파수변화
- 그 외 전원 교란

전원 문제가 의심되는 경우 전원 품질 감사를 받아볼 것을 권장합니다. 자세한 내용은 해당하는 전력 공급 기관 또는 당사에 문의하십시오.

전기 서비스 사양

다음 표는 전기 서비스에 대한 사양입니다.이 요건에 관하여 궁금한 점이 있는 경우 당사의 해 당 서비스 담당자에게 문의하십시오.

요건	사양
입력 전류	최대 1.5A
입력 전압	100 ~ 240 VAC

Nicolet Summit OA 분광기 시작하기

요건	사양
라인 주파 수	50~60
전원 교란	강하,서지 또는 그 외 전원 교란은 입력 전압의 10%를 초과하면 안 됩니 다.
잡음	2 V 미만(일반 모드) 20 V 미만(정상 모드)

Summit 분광기와 개인적 안전

Nicolet Summit OA 분광기는 안전한 장비로 설계되었지만, 정상 사용 시와 유지보수 시 발생할 수 있는 잠재적인 위험으로부터 자신을 보호하려면 몇 가지 예방 조치를 취해야 합니다.

주의 이 설명서에서는 주의



이 설명서에서는 주의해야 할 잠재적인 위험을 소개하지만,이것 이 포괄적인 안내는 아닙니다.이 장비를 사용하기 전에,분광기의 작업장 및 안전 정보를 읽고 이러한 잠재적인 위험에 대해 전반적 으로 이해하십시오.

정상 사용 시의 잠재적인 위험

정상 사용 시 대부분의 위험은 다음과 같은 원인으로 인해 발생합니다.

- 잠재적으로 위험한 샘플과 용제
- 셀렌화 아연(ZnSe)에 대한 노출
- 뜨겁거나 열을 방출하는 분광기의 부품
- 장비의 레이저

잠재적인 위험을 이해하고 필요한 예방 조치를 취하면 가능한 위험이나 장비의 손상을 피할 수 있습니다.

위험한 샘플과 용제

가압 가스나 부식성 또는 가연성 용제 등 잠재적으로 위험한 샘플이나 용제를 사용하거나 측정 하려는 경우 특별한 주의를 기울이십시오.

적절한 환기 사용

분광기에 대한 특별한 환기 요건은 없지만, 특정 유형의 분석 시 추가적으로 환기하는 것이 필 요할 수 있습니다. 독성이 높은 샘플을 분석하거나 적외선 소스와 접촉할 수 있는 용제에 샘플 을 용해하거나 가연성 가스를 샘플링하는 경우 적절한 환기가 되도록 하십시오.

할로겐화 탄화수소가 함유되어 있는 용제를 열분해하면 염산(HCI), 불화 수소산(HF) 또는 포스 겐(COCl₂)이 생성될 수 있습니다.

경고



유독성 물질 흡입 금지. 염산, 불화 수소산, 포스겐은 매우 강한 독 성 물질입니다. 할로겐화 탄화수소를 함유하고 있는 용제를 사용 하는 경우 작업 구역을 적절하게 환기하십시오.

휘발성/가연성 용제

분광기 안의 적외선 소스는 가연성/휘발성 샘플과 용제를 발화시킬 수 있습니다.가연성 샘플 과 용제를 사용하는 경우 다음과 같은 조치를 취하십시오.

- sample compartment window 을 설치한 상태로 작업하십시오.
- 작업 공간에 불꽃이나 다른 점화원이 없고 장비 주변의 대기에 가연성 증기가 모이지 않도 록 방지하는 활성 환기 시스템을 갖추어 적절한 환기가 되도록 하십시오.
- 가연성 용제나 샘플을 장비 근처에 두지 마십시오.
- 가연성 용제나 샘플을 sample compartment에 필요한 시간 이상으로 두지 마십시오.
- 깨끗하고 건조한 공기나 질소로 분광기를 퍼지하십시오.

부식성 용제

sample compartment에서 HCI 또는 HF 증기를 생성하는 용제는 시스템을 심하게 손상시킬 수 있습니다. 할로겐화 용제를 사용하는 경우 깨끗하고 건조한 공기 또는 질소로 장비를 퍼지하십시 오.

알림

장비를 퍼지하지 못하여 발생하는 장비 손상에는 보증이 적용되지 않습니다.

생물학적 위험 또는 방사성 물질과 감염원

조직,체액,감염원,혈액 등의 생물학적 샘플은 전염병을 옮길 수 있습니다.감염성이 있는 물질 을 사용할 때 귀사의 생물학적 안전 프로그램 프로토콜을 따르십시오.

유독성 물질

Nicolet Summit OA 분광계에는 ZnSe(셀렌화 아연) ATR Crystal이 포함되어 있으며, 다른 Summit 장비는 옵션인 ZnSe sample compartment window를 사용할 수 있습니다.

Thermo Scientific

경고



유독성 물질 흡입 금지.

셀렌화 아연(ZnSe)은 유독성입니다. 취급과 노출 관리에 대해서는 www.specac.com에서 제조업체의 ZnSe 안전보건자료를 참조하십시오.

열 원

정상 사용 시 이 분광계의 부품이 상당히 뜨거워질 수 있습니다.분광계의 적외선원과 환기구 주변에 주의하십시오.

바깥쪽을 향하는 적외선원의 표면은 장비 하단에 있으며, 매우 뜨거워질 수 있습니다. 사용 시 또는 사용 직후에는 장비 바닥에 접촉하지 마십시오.

분광기의 환기구는 장비의 왼쪽에 있습니다.정상 사용 시 분광기는 내부 구성부품에서 분광기 바깥으로 이러한 환기구를 통해 뜨거운 공기를 방출합니다.뜨거운 공기가 방출되도록 장비 환 기구 주위에 충분한 공간을 두십시오.

레이저및광학안전

분광기를 정상적으로 사용하고 있을 때는 위험한 수준의 레이저 방사선에 노출되는 일이 없습 니다.서비스 절차 중에 덮개를 분리한 경우 보호복 착용 등의 특별 예방 조치를 취해야 합니다. 필요한 경우 서비스 담당자가 이를 알려 줍니다.

경고



사람 부상 방지.

레이저 빔이나 그 반사물을 응시하지 마십시오.결함 있는 레이저를 교환하 는 경우에도 절대 레이저를 건드리지 마십시오.그러면 레이저 광선 또는 고 압에 노출될 수 있습니다.

유지보수 시 잠재적인 위험

장비 유지보수 작업 시에는 정상 사용 때와 다른 위험에 노출될 수 있습니다.유지보수 시 주된 위험에는 장비 퍼지,장비 내부 구성부품 작업 등이 포함됩니다.

장비 퍼지

특히 습한 환경에서는 깨끗하고 건조한 공기나 질소 공급 장치를 설치하여 분광기를 퍼지하는 것이 좋습니다.장비를 퍼지하면 습한 환경이나 부식성 용제가 유발하는 손상으로부터 내부 광 학 구성부품을 보호할 수 있으며,더 정확한 결과도 확보할 수 있습니다. 분광기의 퍼지 킷을 구매하고 설치하는 자세한 방법은 <u>"퍼지 킷 설치 및 유지보수"</u> 항목을 참조 하십시오.

위험	
	화재 또는 폭발 위험 방지.
\wedge	• 건조한 공기나 질소만 사용하여 분광기를 퍼지하십시오.
	 가연성, 연소성 또는 유독성 가스를 사용하여 이 장비를 퍼지하지 마십 시오. 퍼지 가스에는 오일과 기타 반응성 물질이 없어야 합니다. 소스에 서나 혹은 레이저 흡수로 발생하는 열이 퍼지 가스에서 가연성 가스 또 는 반응성 물질을 발화시킬 수 있습니다.

내부 구성부품 작업

일반적으로, 사용자가 장비 덮개를 분리하거나 내부 구성부품을 교환할 이유는 없습니다. 그러 나 유지보수를 위해 덮개를 분리해야 하는 경우 감전되거나 화상을 입거나 레이저 광선에 노출 될 수 있다는 점을 유념하십시오.



건조제 교환

건조제실을 여는 경우 가연성 액체 또는 가스가 건조제실에 들어오지 않게 막아야 합니다.건조 제 교환 지침은 <u>"건조제 교환"</u> 항목을 참조하십시오.



IR 소스 교환

사용 시 IR 소스는 극도로 뜨거워집니다. IR 소스를 교환해야 하는 경우 화상과 폭발 위험을 방 지하십시오.



분광기에서 소스를 제거하기 전에,장비 전원을 끄고,전원 코드를 분리하 고,모든 퍼지 라인을 분리하고,시스템에서 샘플을 제거하십시오.가연성 액체 또는 가스가 소스실에 들어가면 폭발이 발생할 수 있습니다.이러한 경 우 즉시 당사에 연락하고,이 조건이 해소될 때까지 장비의 전원을 켜지 마 십시오.

요약

Nicolet Summit OA 분광기는 안전하고 견고한 장비이지만, 사용하거나 유지보수할 때 위험에 노 출될 수 있습니다. 정상 사용 시, 잠재적으로 위험한 샘플과 용제를 취급할 때 주의를 기울이고 뜨겁거나 열을 방출하는 분광기의 부품에 접촉하지 말아야 합니다. 유지보수 시, 장비를 퍼지 하거나 내부 구성부품을 취급하거나 건조제를 교환할 때 발생할 수 있는 손상 또는 위험을 피하 려면 예방 조치를 취하십시오.

분광기 포장 풀기 및 시작하기

Thermo Scientific Nicolet Summit OA 분광기의 포장을 풀고 설정하려면 작업 공간에 장비를 놓고, 분광기를 연결하여 전원을 켜고, 간단한 성능 테스트를 실행해야 합니다.

시작하기 전에

알 림

플라스틱 제품 백에서 분광기를 꺼내기 전에 내부 광학장치를 손상시킬 수 있는 응축이 형성 되지 않도록 24시간 동안 기다리십시오.장비가 실온에 도달하기 전에 밀봉된 플라스틱 백을 개봉하여 발생한 손상에는 보증이 적용되지 않습니다.

주의: 분광기의 작업장 및 안전 정보를 검토하여 작업 공간이 장비에 안전하고 적합한지 확인하 십시오.

포장 풀기 및 설치

분광기 포장 풀기/설치 단계는 구매한 옵션에 따라 다소 다릅니다.

1. 제품 상자의 포장을 푸십시오.

제품 상자와 백에서 분광기를 꺼내고 추가 포장을 푸십시오.

이 분광기는 ATR 크리스탈 위에 탈착식 먼지 덮개가 장착되어 제공됩니다.장비를 보관하 거나 이동할 때에는 먼지 덮개를 사용하고 정기적으로 사용할 경우에는 탈거하십시오.

- 2. 전원공급장치를 분광기와 적절한 접지가 되어 있는 AC 전원에 연결하십시오.
- 3. 모니터, 키보드, 마우스를 연결하십시오.
 - a. 모니터를 분광기의 Mini DisplayPort 연결부에 연결하십시오.
 - b. 분광기의 USB 포트 중 하나에 키보드와 마우스를 연결하십시오.

참고 일부 USB와 Mini DisplayPort 기기는 무선 키보드와 마우스 등의 무선 USB 기기에서 나오는 신호를 방해하는 것으로 알려져 있습니다. 무선 키보드와 마우스를 사용 할 때 입력 지연 또는 누락 등의 성능 문제가 발생하는 경우 제공된 USB 연장 케이블 을 사용하여 무선 USB 수신기를 USB 포트에서 멀리 두십시오.

4. 장비의 전원 버튼을 눌러서 분광기를 켜십시오.장비 예열 중에 분광기의 LightBar에 순환되

는 녹색 라이트가 켜지고,장비 사용 준비가 완료되면 고정된 녹색 막대가 표시됩니다.

장비를 켜거나 끌 때 내부 폴리스티렌 기준 물질이 움직이는 소리가 들릴 수 있습니다.이것 은 정상입니다.



분광기를 처음 켜면 새 암호를 입력하라는 프롬프트 메시지가 표시됩니다.이 때는 암호를 입력 하지 않는 것이 좋습니다.암호를 입력하지 않고 진행한 후 설정을 완료하면 조직의 정책에 따 라 장치의 계정과 암호를 관리하십시오.

참고 시스템을 켰을 때 화면이 검은색이면 모니터 전원, DisplayPort 연결, 모니터 입력을 확 인하십시오. 입력 설정 변경에 대한 자세한 내용은 모니터와 함께 제공된 설명서를 참조하 십시오.

5. OMNIC Paradigm 소프트웨어를 여십시오.

소프트웨어가 시작되고 장비에 연결하는 데 몇 분 정도 걸릴 수 있습니다.소프트웨어가 화 면의 오른쪽 상단에 연결 상태를 표시합니다.상태에 연결된 것으로 표시되면 다음 단계로 진행하십시오.



참고 장비를 연결하기 전에 소프트웨어에서 메뉴를 열면 Align(조절)과 Calibrate(보정)가 비활성화됩니다.이를 활성화하려면 장비가 연결되어 있을 때 도구 모음의 Dashboard(대 시보드)를 클릭하십시오.

6. 분광기의 성능을 확인하십시오.

- a. 분광기 광학장치를 조절하십시오.
 - i. Desktop(데스크탑)인터페이스에서 Acquire Data(데이터 획득) > Diagnostics(진단) > Align Spectrometer(분광기 조절)로 이동하십시오.
 - ii. Start(시작)를 선택하여 조절을 시작하십시오.조절이 정상적으로 완료되면 화면에 'Align complete(조절 완료)'메시지가 표시됩니다.
- b. 레이저를 보정하십시오
 - i. Acquire Data(데이터 획득) > Diagnostics(진단) > Laser Calibration(레이저 보정)으로 이동하십시오.
 - ii. Start(시작)를 선택하여 보정을 시작하십시오.
 - iii. 레이저 보정이 완료될 때까지 몇 분 정도 걸릴 수 있습니다. 보정이 완료되면 디스플 레이에 'Measurement complete(측정 완료)' 메시지가 표시되고, 레이저 주파수의 변화 가 표시됩니다.

장비가 열평형에 도달할 때까지 최대 12시간 정도 걸립니다.최고의 결과를 얻으려면 처음 12시간 작동 후 조절과 보정을 다시 실행하십시오.

c. Nicolet Summit OA - PV Test(Nicolet Summit OA - PV 테스트) 작업 플로우를 실행하십시오.

성능 확인 작업 플로우에서는 기준 물질을 사용하여 일련의 표준 테스트를 실행하고 장 비 작동을 확인하고 데이터의 정확성을 확보합니다.

- i. Desktop(데스크탑)인터페이스의 Dashboard(대시보드)에서 Workflows(작업 플로우) 창으로 스크롤 이동하십시오.
- ii. Workflow(작업 플로우) 창에서 Nicolet Summit OA PV Workflow(Nicolet Summit OA PV 작업 플로우)를 마우스 오른쪽 버튼으로 클릭하고 Run(실행)을 선택하십시오.

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

- iii. 화면에 표시되는 메시지에 따라 작업 플로우를 완료하십시오.
- iv. PV Results(PV 결과) 표가 표시되면 Finish(마침)를 선택하여 작업 플로우를 완료하고 대시보드로 돌아가거나, Measure Squalane Spectrum(스쿠알란 스펙트럼 측정)을 선 택하여 작업 플로우의 선택적 스쿠알란 부분으로 진행하십시오. 스쿠알란 스펙트럼 을 측정하면 시스템 성능을 확인하기 위한 추가적인 기준 확인이 제공됩니다.



다음 단계

축하합니다! 장비 사용 준비가 완료되었습니다. 그러나 데이터 수집을 시작하기 전에, 나중에 발생할 수 있는 잠재적인 문제로부터 시스템을 보호할 수 있도록 시스템을 백업하고 복구 드라 이브를 만드십시오. Windows 10에는 데이터를 백업하고 복구할 수 있는 여러 가지 기본 제공 도 구가 들어 있습니다.

복구 드라이브와 시스템 이미지를 만들고 시스템을 복원하는 지침은 <u>"데이터 백업 및 복구"</u> 항 목을 참조하십시오.

분광기로 샘플을 식별하고 분석하는 자세한 방법은 thermofisher.com/ftir-help에서 OMNIC Paradigm 설명서와 튜토리얼을 참조하십시오.

Operator(작업자)인터페이스로 전환하려면 **View / Display(보기/표시) > Operator(작업자)**로 이 동하십시오.

Package Editor(패키지 편집기)와 Operator(작업자) 인터페이스를 사용하여 맞춤 솔루션을 만들고 실행하는 자세한 지침은 OMNIC Paradigm 소프트웨어 사용자 설명서의 'Custom Solutions(맞춤 솔루션)'를 참조하십시오.

데이터 백업 및 복구

데이터 수집을 시작하기 전에, 나중에 발생할 수 있는 잠재적인 문제로부터 시스템을 보호할 수 있도록 사전 주의 단계를 수행해야 합니다. Windows 10에는 데이터를 백업하고 복구할 수 있는 여러 가지 기본 제공 도구가 들어 있습니다.

중요! 하드 드라이브 또는 운영 체제가 손상된 경우 데이터를 복구할 수 있도록 복구 미디어를 만들고, 시스템 이미지를 만들고, 데이터 백업 절차를 설정하십시오.

복구 드라이브 만들기

설치된 Windows가 손상되는 경우에도 시스템 이미지 복구,시스템 복원,시작 설정 등의 Windows 10 고급 시작 옵션에 액세스할 수 있도록 복구 드라이브를 만드십시오.

필요한 품목

• 저장 공간이 8 GB이상인 FAT32 포맷 플래시 드라이브

알림

드라이브의 모든 내용이 이 프로세스에서 삭제됩니다.진행하기 전에, 보관할 모든 파일을 백 업하십시오.

◆ 복구 드라이브를 만들려면

- 1. 분광기 측면의 USB 포트 중 하나에 플래시 드라이브를 삽입하십시오.
- Windows 10 작업 표시줄의 검색 상자에서 'Create Recovery Drive'를 검색하고 Create Recovery Drive(복구 드라이브 만들기)를 선택하십시오.



3. 장치 변경을 허용할 것인지 물으면 Yes(예)를 선택하십시오.

- 4. 'Back up system files to the recovery drive(복구 드라이브에 시스템 파일 백업)'의 선택을 지우고 Next(다음)를 클릭하십시오.
- 5. 가용 드라이브 목록에서 원하는 드라이브를 선택하고 Next(다음)를 클릭하십시오.

 	×
Select the USB flash drive	
The drive must be able to hold at least 512 MB, and everything on the drive will be deleted.	
Available drive(s) └─ D∿ (RECOVERY)	
Next	Cancel

- 6. Create(만들기)를 클릭하십시오.이 프로세스가 완료되면 Finish(마침)를 클릭하여 창을 닫으십시오.
- 7. 플래시 드라이브를 안전하게 꺼내서 안전한 곳에 보관하십시오.드라이브에 'Summit 복구 드라이브'라고 써 놓으면 좋습니다.

복구 미디어 만들기에 대한 자세한 내용은 Microsoft 온라인 도움말을 참조하십시오.

시스템 이미지 만들기

시스템 이미지는 데이터와 파일,설치된 응용프로그램,기타 설정을 포함한 전체 시스템의 스 냅샷입니다. 전체 하드 드라이브 또는 운영 체제에 장애가 발생한 경우 이 시스템 이미지를 사 용하여 장치를 이미지가 만들어진 때의 상태로 복원할 수 있습니다. 시스템 이미지를 얼마나 자 주 업데이트할 것인가는 사용자의 판단이지만,장비를 처음 설치했을 때의 이미지를 보관하는 것이 좋습니다.

시스템 이미지에 복사하는 데이터의 양에 따라, 프로세스에 시간이 다소 걸릴 수 있습니다. 대 용량의 데이터를 복사하는 경우 프로세스가 완료될 때까지 1시간 이상이 걸릴 수 있습니다.

필요한 품목

- NTFS 포맷 SSD
- ◆ 시스템이미지를 만들려면
- 1. 분광기의 USB 포트 중 하나에 SSD를 연결하십시오.
- 2. Windows 10에서 제어판을 열고 백업 및 복원(Windows 7)을 선택하십시오.

3. 왼쪽 패널에서 시스템 이미지 만들기를 선택하십시오.

← → × ↑ 🐌 > Control P	anel > 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. 하드 디스크를 선택하고 목록에서 해당 SSD를 선택하십시오. 다음을 클릭하십시오.

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

- 5. 백업 설정을 확인하고 Start backup(백업 시작)을 클릭하십시오.
- 6. '시스템 복구 디스크를 만드시겠습니까?'라고 물으면 No(아니요)를 클릭하십시오.
- 7. Close(닫기)를 클릭하고 Summit에서 SSD를 안전하게 꺼내십시오. 드라이브에 Summit 복구 라고 써서 안전한 곳에 보관하십시오.

시스템 복원

운영 체제에 장애가 발생한 경우 복구 드라이브와 시스템 이미지를 사용하여 시스템을 복원하 십시오.

필요한 품목

- USB 복구 드라이브
- 시스템 이미지가 들어 있는 SSD
- ◆ 복구 드라이브와 시스템 이미지를 사용하여 시스템을 복원하려면
- 1. Summit 분광기를 끄십시오.
- 2. Summit 분광기의 USB 포트 중 하나에 외부 키보드를 연결하십시오.
- 3. USB 포트 중 하나에 USB 복구 드라이브를 연결하십시오.
- 4. 시스템을 켜십시오.
- 5. 화면에 Thermo Scientific 시작 화면이 표시되면 키보드의 'Esc'를 여러 번 눌러서 시스템 BIOS를 여십시오.

팝업 대화 상자에 '저장하지 않고 종료하시겠습니까?' 메시지가 표시되면 화살표 키를 사용 하여 아니요를 선택하십시오.

- 6. 화살표 키를 사용하여 Save & Exit(저장 후 종료) 탭으로 이동하십시오.
- 7. Boot Override(부팅 다시 정의) 섹션으로 이동하여 해당 복구 플래시 드라이브를 선택하십시 오. Enter를 눌러서 복구 이미지로 부팅하십시오.



- 8. '키보드 레이아웃 선택' 화면이 표시되면 USB 포트에서 복구 플래시 드라이브를 꺼내고 시 스템 이미지 SSD를 삽입하십시오.
- 9. 원하는 키보드 레이아웃을 선택하십시오.
- 10. 옵션 선택 화면에서 Troubleshoot(문제 해결)을 선택하십시오.

Choose an option	
Continue Exit and continue to Windows 10	
Troubleshoot Reset your PC or see advanced options	
Turn off your PC	

- 11. Advanced options(고급 옵션)을 선택하십시오.
- 12. 고급 옵션에서 System Image Recovery(시스템 이미지 복구)를 선택하십시오.

	System Restore		Startup Repair
	Use a restore point recorded on your PC to restore Windows	< \$ >	Fix problems that keep Windows fro loading
	Go back to the		Command Prompt
.	previous version	C:\	advanced troubleshooting
	System Image Recovery		Startup Settings Change Windows startup behavior
	Recover Windows using a specific system image file	- 1	

- 13. Windows 10을 대상 운영 체제로 선택하십시오.
- 14. '시스템 이미지 백업을 선택합니다.'창에서 최신 가용 시스템 이미지를 사용하거나 다른 시 스템 이미지를 선택하십시오. Next(다음)을 클릭하십시오.

🍋 Re-image your computer			×
	Select a syst This computer will b Everything on this information in the s	em image backup e restored using the system image. computer will be replaced with the system image.	
	Troubleshooting inf http://go.microsoff Ouse the latest a	formation 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	
		< Back Next > Can	cel

Thermo Scientific

15. Next(다음)을 클릭하여 추가적인 복원 옵션을 기본 상태로 두고, 마침을 클릭하여 시스템 이 미지를 복원하십시오. 예를 클릭하여 확인하십시오.

복원에는 약 15분 정도가 소요되며, 완료되면 시스템이 자동으로 다시 부팅됩니다.

시스템이 다시 켜지면 분광기에서 SSD를 안전하게 꺼내고 복구 드라이브와 시스템 이미지 SSD를 안전한 곳에 보관하십시오.

시스템을 이미지로 복원하는 방법은 Microsoft 온라인 도움말을 참조하십시오.

시스템 데이터 백업

복구 미디어와 시스템 이미지를 만드는 것 외에,정기적으로 데이터를 백업해야 합니다. Windows 10에는 파일 히스토리 도구를 포함하여 파일을 백업하는 여러 가지 옵션이 있습니다. 사용자 조직의 정보 기술(IT) 담당자는 조직의 필요에 가장 적합한 백업 방법이 무엇인지 결정 해야 합니다.

USB 무선 어댑터 설치

옵션 품목인 USB 무선 어댑터를 사용하여 Nicolet Summit OA 분광기를 무선 네트워크에 연결할 수 있습니다.분광기와 함께 어댑터를 구매한 경우 미리 설치되어 제공됩니다.

USB 무선 어댑터를 설치하려면 어댑터를 아무 USB 포트에 삽입하십시오. 영구적으로 설치하 려면 장비 덮개 아래의 포트에 어댑터를 설치하십시오.

◆ USB 무선 어댑터를 설치하려면

1. 어댑터를 빈 USB 포트에 삽입하십시오.



2. 시스템이 자동으로 어댑터를 인식합니다.어댑터를 사용하여 네트워크에 연결하는 방법은 Microsoft 지원 자료를 참조하십시오.

◆ USB 포트를 사용하여 장비 덮개 아래에 설치하려면

이렇게 설치하려면 크리스탈 플레이트와 장비 덮개를 탈거해야 합니다.

- 1. "덮개 교환"의 지침에 따라 크리스탈 플레이트와 장비 덮개를 탈거하십시오.
- 2. 덮개를 탈거한 상태에서 어댑터를 빈 USB 포트에 삽입하십시오.

Nicolet Summit OA 분광기 시작하기



3. 장비 덮개와 크리스탈 플레이트를 다시 장착하십시오.

이제 장비의 무선 네트워크 연결 준비가 완료되었습니다.

분광계 LightBar 이해

Thermo Scientific Nicolet Summit OA 분광기에는 제품 품질과 장비 상태를 한눈에 볼 수 있는 신속 하고 시각적인 통과/실패 표시를 제공하는 편리한 LightBar가 있습니다.

다음 표는 모든 LightBar 신호를 보여 줍니다.

예 열



기존 수집

신호	시스템 상태	설명
고정된 녹색	사용 준비 완료	전체 LightBar가 고정된 녹색을 표시합니다.
녹색 코멧	데이터 수집 중	녹색 라이트의 작은 부분이 앞뒤 로 움직입니다.

스마트 수집

신호	시스템 상태	설명
파란색 펄스	스마트 배경 수 집 중	LightBar 전체가 파란색으로 변합 니다.
파란색 코멧	샘플 데이터 수 집 중(사용자가 시작함)	파란색 라이트의 작은 부분이 앞 뒤로 움직입니다.

분석 결과

신호	시스템 상태	설명
녹색 % 채움	일 치 값 또는 임 계 값을 초과 하 는 QCheck 결 과	일치 값을 반영하도록 녹색이 채 워집니다. 예를 들어, 일치 값이 90이라면 약 90%를 채웁니다.
오렌지색 % 채움	일 치 값 또 는 임 계 값 미 만 인 QCheck 결 과	일치 값을 반영하도록 오렌지색 이 채워집니다.예를 들어,일치 값이 30 이라면 약 30% 를 채웁니 다.

진 단

신호	시스템 상태	설명
빨간색 깜박임	시스템 오류	LightBar 전체가 빨간색으로 깜박 입니다. 오류에 대한 설명은 OMNIC Paradigm 소프트웨어의 System Status(시스템 상태)를 참조하십시오.

OMNIC Anywhere 앱으로 데이터 보기

OMNIC Anywhere는 연결된 PC, Apple 컴퓨터, Android 또는 iOS 기기에서 데이터를 보거나 탐색 하거나 공유할 수 있는 클라우드 기반 응용프로그램입니다.

2-1 그림:웹 브라우저의 OMNIC Anywhere



Nicolet Summit OA 분광기와 무료 연결 계정을 사용하여 강의실 또는 실험실의 샘플을 측정하고, 연결 계정에 데이터를 업로드하고, 사무실 또는 연구실의 다른 장치에 있는 데이터를 보거 나 탐색하거나 공유할 수 있습니다.

연결 계정을 만들고 OMNIC Anywhere로 데이터를 보는 지침은 '<u>클라우드 지원 FTIR 분광법</u>'을 참조하십시오. [별도 공란]

옵션 및 액세서리

Nicolet Summit OA를 옵션 품목인 퍼지 킷과 함께 사용할 수 있습니다.

퍼지	₹I	설치	및	유지보수	32
	\sim		~	가지포 [

퍼지 킷 설치 및 유지보수

건조한 공기 또는 질소로 분광기를 퍼지하면 습기나 다른 환경 오염으로부터 내부 구성부품을 보호할 수 있습니다.퍼지 킷을 설치하려면 밸브와 조절기를 조립하고,건조제 카트리지를 교 환하고,퍼지 가스를 연결하고,압력과 유속을 설정해야 합니다.

폭발 위험 방지



가연성,연소성 또는 유독성 가스를 사용하여 이 장비를 퍼지하지 마십시오. 퍼지 가스에는 오일과 기타 반응성 물질이 없어야 합니다.소스에서나 혹은 레이저 흡수로 발생하는 열이 퍼지 가스에서 가연성 가스 또는 반응성 물질 을 점화시킬 수 있습니다.건조한 공기나 질소만 사용하여 장비를 퍼지하십 시오.

알 림

경고

항상 씰과 건조를 유지하거나 장비를 퍼지하는 것이 좋습니다.장비를 올바르게 건조하거나 퍼지하지 못해서 발생한 손상에는 보증이 적용되지 않습니다.이 요건에 관하여 궁금한 점이 있는 경우 당사에 문의하십시오.

필요한 설정

퍼지 킷을 설치하기 전에, Nicolet[™] Summit 분광기에 대한 Thermo Scientific 사양에 맞는 건조 공 기원 또는 질소가 필요합니다. 자세한 내용은 장비의 작업장 및 안전 설명서를 참조하십시오.

참고 최선의 결과를 얻으려면 퍼지 가스를 이슬점 -70 °C(-94 °F) 이하로 건조시키십시오.

필요한 품목

퍼지 킷 외에,다음 도구와 품목이 필요합니다.

- 3/4인치 양구 렌치
- 11/16인치 양구 렌치
- #2 십 자 드 라 이 버
- 3/32인치 육각 키
- 7/64인치 육각 키
- 스레드씰 테이프('배관용 테이프' 또는 'Teflon 테이프')

◈ 퍼지 킷을 설치하려면

- 1. 배관 어셈블리를 퍼지 가스 소스에 부착하십시오.
 - a. 조절기 밸브와 1/4인치 수 장착 장치 또는 3/8인치 암 장착 장치를 퍼지 가스 소스에 설치 하십시오. (해당 퍼지 가스 소스에 적합한 밸브와 장착 장치를 선택하십시오.)



b. 1/4인치 수 장착 장치를 사용하는 경우 다음 단계로 진행하십시오.

퍼지 가스 소스에 3/8인치 암 장착 장치를 사용하는 경우 퍼지 킷에 함께 제공된 3/8인 치~1/4인치 축소 니플을 설치하십시오. 설치하기 전에 스레드씰 테이프를 축소 니플에 감고, 11/16인치 양구 렌치를 사용하여 연결부를 조이십시오.



c. 스레드씰 테이프를 축소 니플 또는 1/4인치 수 장착 장치에 감은 후 압력 연결 장치를 설 치하십시오. 3/4인치 양구 렌치를 사용하여 연결부을 조이십시오.



d. 벽면 배관 어셈블리의 수 입구를 빠른 해제 장착 장치에 단단히 끼우십시오.

- 2. 분광기를 끄십시오.
 - a. 분광기를 끄고, 전원 버튼을 누르십시오.
 - b. 전원 코드를 분리하십시오.
- 3. "크리스탈 플레이트 탈거"
- 4. 장비 덮개를 탈거하지 마십시오.
 - a. 장비가 열 핀에 놓이도록 장비를 조심스럽게 뒤로 기울이십시오.
 - b. 7/64 육각 키를 사용하여,덮개를 장비 바닥판에 고정하고 있는 나사 4개를 푸십시오.



- c. 장비를 피트에 조심스럽게 뒤로 기울이십시오.기기를 뒤로 젖히면 4개 나사가 빠집니다. 잃어버리지 않도록 주의하십시오.
- d. 덮개를 바닥판에서 똑바로 들어 올리십시오.



5. 건조제 카트리지를 분리하십시오.

a. #2 십자 드라이버를 사용하여 건조제 카트리지에서 두 계류 나사를 풀고, 카트리지를 수 직으로 들어 올려서 장비에서 꺼내십시오.



b. 큰 O 링을 분리하여 버리십시오.



참고 O링을 건조제실에 떨어뜨리지 않도록 주의하십시오.

- 6. 퍼지 킷과 밀봉된 건조제 패킷을 여십시오.
- 7. 패킷에서 O 링을 분리하고 장비 바닥판에 설치하십시오(이전 그림 참조). 새 O 링을 제자리 에 눌러서(굴리지 않음) 홈에 올바르게 장착되도록 하십시오.
- 8. 패킷에서 건조제 카트리지와 퍼지 장착 장치를 분리하고, 건조제실에 카트리지를 삽입하

고, 0 링 위로 올바르게 장착되었는지 확인한 후, 두 나사를 사용하여 고정하십시오.

카트리지의 올바른 방향을 기억하십시오.카트리지가 설치되면 장비 앞쪽에서 카트리지의 레이블을 읽을 수 있습니다.



- 9. 퍼지 라인을 기기 뒷면의 퍼지 입구로 통과시키십시오.
 - a. 퍼지 튜브를 덮개 슬롯으로 라우팅하면서 덮개를 받침대 위로 조심스럽게 내리십시오.



- b. 덮개를 완전히 내리십시오.덮개를 제자리로 내리면 퍼지 튜브가 덮개와 방열판 사이의 슬롯에 끼이게 됩니다.
- c. 퍼지 튜브를 옆으로 밀고 장비를 뒤쪽으로 기울여 열 핀에 올려 놓으십시오.
- d. 이전에 분리한 4개 나사를 삽입하십시오.모든 나사를 조이십시오.
- e. 장비를 피트에 뒤로 기울이십시오.
- f. 남은 길이의 튜브는 장비에 밀어 넣거나 필요에 따라 빼내십시오.



- 10. "크리스탈 플레이트 교환".
- 11. flow coupler를 벽면 배관 어셈블리('분광기 연결'표시)에서 퍼지 튜브로 연결하십시오.
- 12. 퍼지 가스 컨트롤을 설정하십시오.
 - a. 주 조절기 밸브를 열고 아래와 같이 퍼지 가스 컨트롤을 설정하십시오(자세한 내용은 퍼지 가스 컨트롤 설정 참조).

하드웨어	압 력 (psig)	유속(scfh)
Summit 분광기	5	1

b. 전원 코드를 장비에 연결하고 전원을 켜십시오.

c. 장비가 완전히 퍼지되도록 30~60분 정도 기다리십시오.

퍼지 가스 컨트롤 설정

퍼지 가스 컨트롤을 올바르게 설정하면 진동을 유발하지 않고 습도로부터 분광기를 보호할 수 있습니다.

최선의 결과를 얻으려면 퍼지 가스를 이슬점 -70°C(-94°F) 이하로 건조시켜야 합니다.

◆ 퍼지 가스 컨트롤을 설정하려면

- 1. 주 밸브를 열어서 조절기를 통하는 퍼지 가스 흐름을 시작하십시오.
- 2. 게이지의 압력이 5 psig(34 kPa)가될 때까지 분광기의 압력 조절기를 조절하십시오.
- 3. 분광기의 유량계를 1 scfh(0.47 l/분)로 설정하십시오.

알림

A 유속이 권장 값을 초과하면 진동이 발생하여 데이터 품질에 영향을 미칠 수 있습니다.

퍼지 가스 필터 확인 및 교환

퍼지 필터가 노란색이거나 다른 색으로 변색되거나 잔류물 또는 이물질로 오염된 경우 필터를 교환하십시오.

알 림

항상 씰과 건조를 유지하거나 장비를 퍼지하는 것이 좋습니다. 씰과 건조를 유지하거나 퍼징 하지 못한 결과로 인한 장비 손상에는 보증이 적용되지 않습니다. 이 요건에 관하여 궁금한 점이 있는 경우 당사에 문의하십시오.

3-1 그림: 퍼지 필터는 압력 게이지 아래의 플라스틱 보울 안에 있습니다.



참고 부품을 주문하려면 기술 지원팀에 문의하십시오.

◆ 퍼지 가스 필터를 교환하려면

- 1. 주 밸브의 퍼지 가스를 끄십시오. 유량계 또는 압력 조절기를 낮추지 마십시오.
- 2. 필터가 들어 있는 플라스틱 보울을 분리한 후 필터를 분리하십시오. (둘 모두 손으로 풀 수 있습니다.)



- 3. 새 필터를 설치한 후 보울을 다시 설치하십시오.
- 4. 주 밸브를 열어서 장비로 퍼지 흐름을 켜고 퍼지 가스 유속이 적절한지 확인하십시오.

4 장

유지보수

Nicolet Summit OA FTIR 분광기는 관리가 거의 필요 없으며, 일일 유지보수는 전혀 필요 없습니 다. 그러나 기본적인 관리를 해 주면 지속적으로 최고 성능을 유지할 수 있습니다.

Summit 분광기의 유지보수 일정	42
분광기 광학장치 정렬	44
레이저 주파수 보정	45
장비와 크리스탈 청소	46
크리스탈 플레이트 탈거 및 교환	48
덮개 교환	
건조제 교환	54
소스 교환	60

Summit 분광기의 유지보수 일정

Thermo Scientific Nicolet Summit OA FTIR 분광기는 관리가 거의 필요 없으며, 일일 유지보수는 전혀 필요 없습니다.그러나 기본적인 관리를 해 주면 지속적으로 최고 성능을 유지할 수 있습 니다. 이 항목의 지침을 따르십시오. 최고의 성능을 유지하려면 분광기의 전원을 켜 두십시오.

알림

정전기는 분광기의 중요 구성부품을 영구적으로 손상시킬 수 있습니다.이러한 손상을 방지 하려면 다음 권장 사항을 따르십시오.

- 전원공급장치를 분리하기 전에, 분광기의 금속 베이스를 접촉하여 축적되어 있을 수 있는 정전기를 모두 방전하십시오.
- 교환 부품은 장비에 설치할 준비가 완료될 때까지 포장 상태 그대로 두십시오.

주간 유지보수

SPECTROMETER 성능 확인

OMNIC Paradigm 소프트웨어에는 분광기에 대한 성능 확인 작업 플로우가 포함됩니다. PV 작업 플로우는 일련의 표준 테스트를 실행하여,장비 작동을 확인하고 데이터의 정확성을 확보합니 다.

주 1회 이상, PV 작업 플로우를 실행하는 것이 좋습니다.

분광기 청소

이 분광기는 권장된 대로만 청소할 수 있습니다. "분광기 청소" 항목을 참조하십시오.

월간 유지보수

습도 표시기 확인

이 분광기의 광학 구성부품은 공기 중의 과도한 습기에 의해 쉽게 손상될 수 있습니다.이 분광 기는 밀봉되어 있으며,그 구성부품은 습기를 흡수하는 두 건조제 캐니스터에 의해 보호됩니 다. OMNIC Paradigm 소프트웨어에서 장비 습도를 모니터링하십시오.

알 림

항상 분광기를 밀봉하고 건조한 상태로 유지하거나 밀봉하고 마른 공기 또는 질소로 퍼지하 는 것을 권장합니다. 씰과 건조를 유지하거나 퍼징하지 못한 결과로 인한 장비 손상에는 보증 이 적용되지 않습니다. 이 요건에 관하여 궁금한 점이 있는 경우 당사에 문의하십시오.

퍼지 가스 필터 확인

분광기를 질소나 건조한 공기로 퍼지하는 경우 월 1회 이상 퍼지 필터를 확인하십시오.

필터가 노란색이거나 다른 색으로 변색되거나 잔류물 또는 이물질로 오염된 경우 필터를 교환 하십시오.자세한 내용은 "퍼지 가스 필터 확인 및 교환" 항목을 참조하십시오.

4-1 그림: 퍼지 필터는 압력 게이지 아래의 플라스틱 보울 안에 있습니다.



분광기 광학장치 정렬

Nicolet Summit OA - PV Test(Nicolet Summit OA - PV 테스트) 작업 플로우에서 테스트 실패가 표 시되는 경우 분광기 광학장치를 조절하십시오. 분광기를 조절하면 검출기에 도달하는 에너지 를 최적화하고 검출기 신호를 극대화할 수 있습니다. 성능 테스트에 실패하면 분광기의 광학장 치를 조절하고 레이저 주파수를 보정하고 다시 테스트를 실행하십시오. 여전히 테스트에 실패 하면 현지 기술 지원 담당자에게 도움을 요청하십시오.

◆ 분광기 광학장치를 조절하려면(터치스크린 인터페이스)

- 1. 홈 화면에서 Diagnostics(진단) 아이콘())을 선택하여 Diagnostics(진단) 보기를 여십시오.
- 2. Align(조절) 탭을 열고 Align(조절)을 선택하십시오.

시스템이 자동으로 분광기를 조절하고 조절이 완료되면 이를 표시합니다.

♦ 분광기 광학장치를 조절하려면(작업자 인터페이스)

- 1. Administrator(관리자) 영역으로 이동하고 Diagnostics(진단) 아이콘(⁽⁾)을 선택하여 Diagnostics(진단) 보기를 여십시오.
- 2. Align(조절) 탭을 열고 Align(조절)을 선택하십시오.

시스템이 자동으로 분광기를 조절하고 조절이 완료되면 이를 표시합니다.

- ♦ 분광기 광학장치를 조절하려면(데스크탑 인터페이스)
- 1. Acquire Data(데이터 획득) > Diagnostics(진단) > Align Spectrometer(분광기 조절)를 선택하 십시오.
- 2. Align Spectrometer(분광기 조절) 대화 상자에서 Start(시작)를 클릭하십시오.

시스템이 자동으로 분광기를 조절하고 조절이 완료되면 이를 표시합니다.

레이저 주파수 보정

Nicolet Summit OA - PV Test(Nicolet Summit OA - PV 테스트) 작업 플로우에서 테스트 실패가 표 시되는 경우 레이저 주파수를 보정하십시오. 테스트에 실패하면 먼저 분광기를 조절한 후 레이 저 주파수를 보정하고 다시 테스트를 실행하십시오. 여전히 테스트에 실패하면 현지 기술 지원 담당자에게 도움을 요청하십시오.

- ♦ 레이저를 보정하려면(터치스크린 인터페이스)
- 1. 홈 화면에서 Diagnostics(진단) 아이콘()을 선택하여 Diagnostics(진단) 보기를 여십시오.
- 2. Laser Calibration(레이저 보정) 탭을 열고 Calibrate(보정)를 선택하여 보정을 시작하십시오.

레이저 보정이 완료되면 레이저 주파수가 변경되었는지 여부를 알려 주는 메시지가 표시되고, 새 주파수가 표시됩니다.

- ♦ 레이저를 보정하려면(작업자 인터페이스)
- 1. Administrator(관리자) 영역으로 이동하고 Diagnostics(진단) 아이콘()을 선택하여 Diagnostics(진단) 보기를 여십시오.
- 2. Laser Calibration(레이저 보정) 탭을 열고 Calibrate(보정)를 선택하여 보정을 시작하십시오.

레이저 보정이 완료되면 레이저 주파수가 변경되었는지 여부를 알려 주는 메시지가 표시되고, 새 주파수가 표시됩니다.

- ◆ 레이저를 보정하려면(데스크탑 인터페이스)
- 1. Acquire Data(데이터 획득) > Diagnostics(진단) > Laser Calibration(레이저 보정)을 선택하 십시오.
- 2. Laser Calibration(레이저 보정) 대화 상자에서 Start(시작)를 클릭하여 레이저를 보정하십시 오.

시스템이 레이저 보정을 완료하면 주파수가 변경되었는지 여부를 알려 주는 메시지가 표시 되고,새 주파수가 표시됩니다.

장비와 크리스탈 청소

Thermo Scientific Nicolet Summit OA 분광기를 청소하기 전에, 분광기를 끄고 전원공급장치를 분 리하십시오.

분광기 청소

부드럽고 깨끗한 천에 연한 용제(예: Windex)를 적셔서 장비 외부를 부드럽게 청소하십시오.천 에 뿌려서 외부의 도장면만 닦으십시오.장비 뒷면의 전자장치에 용제가 묻지 않도록 주의하십 시오.

먼지는 장비 뒤쪽의 전자장치 인클로저에 쌓여서 잠재적으로 열 방출을 방해할 수 있으며,이 때문에 전자 구성부품의 수명이 단축될 수 있습니다.

장비 뒤쪽의 먼지를 제거하려면 압축 공기를 사용하여 먼지를 불어 내십시오. 분광기 뒤쪽의 먼 지를 제거할 때 일체 액체를 사용하지 마십시오.



거친 세제,용제,화학 물질 또는 연마제를 사용하지 마십시오.마감이 손상될 수 있습니다.

ATR 크리스탈 플레이트 청소

깨끗하고 마른 비 마모성 천이나 티슈를 사용하여 ATR 크리스탈을 청소하십시오.



- 깨끗한 천이나 티슈를 사용하여 크리스탈에서 액체를 닦아내십시오. 사용한 티슈는 버리십 시오.
- 일회용 피펫을 사용하여 크리스탈과 크리스탈 주변 영역에 헵탄 몇 방울을 떨어뜨리십시 오.
- 깨끗한 새 티슈로 크리스탈의 길이를 따라 닦아서 액체를 닦아내십시오. 크리스탈 골 주변 을 닦으십시오.



4. 필요한 대로 반복하고, 항상 깨끗하고 마른 천이나 티슈로 마무리하십시오.

크리스탈 플레이트 탈거 및 교환

크리스탈 플레이트는 12~18개월 간격으로 또는 필요 시 교환해야 합니다.

유독성 물질 흡입 금지.



셀렌화 아연(ZnSe)은 유독성입니다. 취급과 노출 관리에 대해서는 www.specac.com에서 제조업체의 ZnSe 안전보건자료를 참조하십시오.

크리스탈 플레이트를 청소하거나 다시 장착할 때 보호 장갑을 착용하십시 오.

알 림

경고

덮개가 탈거되었을 때 장비에 물건이 떨어지지 않도록 주의하십시오.

필요한 품목

3/32"육각 키

크리스탈 플레이트 탈거

- 1. 크리스탈에서 먼지 덮개를 탈거하십시오.
- 3/32"육각 키를 사용하여 크리스탈 플레이트를 장비에 고정하고 있는 나사 3개를 완전히 푸 십시오.



- 3. 크리스탈 플레이트를 장비에서 똑바로 들어 올리십시오.
- 깨끗하고 마른 천을 사용하여 플레이트가 놓인 구역 아래 플랜지 주변 영역을 청소하십시 오.

크리스탈 플레이트 교환

- 1. 새 플레이트를 플랜지 위의 제자리에 놓으십시오.
- 2. 와셔를 나사에 끼우고 세 구멍에 나사를 넣으십시오. 나사를 구멍에 똑바로 내려야 합니다.
- 크리스탈 플레이트를 누른 상태에서 플레이트 양쪽에 있는 2개의 나사를 조인 후 장비 뒷면 에 가장 가까운 세 번째 나사를 조이십시오. 손가락으로 모든 나사를 조이십시오.

나사가 자유롭게 회전하는 경우 나사를 위로 올린 후 다시 삽입하여 정렬을 수정하십시오.

덮개 교환

필요한 도구

시작하기 전에,다음 도구를 구비해야 합니다.

- 7/64 육각 키
- 3/32 육각 키

◆ 분광계 덮개를 교환하려면

- 1. 분광기를 끄십시오.
 - a. 분광기를 끄고, 전원 버튼을 누르십시오.
 - b. 전원 코드를 분리하십시오.
 - c. 퍼지를 끄고, 해당하는 경우 장비에 연결된 퍼지 라인을 분리하십시오.
- 2. 크리스탈 플레이트를 탈거하십시오.
 - a. 크리스탈에서 먼지 덮개를 탈거하십시오.
 - b. 크리스탈 플레이트를 장비에 고정하고 있는 나사 3개를 완전히 푸십시오.



Thermo Scientific

- c. 크리스탈 플레이트를 장비에서 똑바로 들어 올리십시오.
- d. 깨끗하고 마른 천을 사용하여 플레이트가 놓인 구역 아래 플랜지 주변 영역을 청소하십 시오.
- 3. 장비 덮개를 탈거하지 마십시오.
 - a. 장비가 열 핀에 놓이도록 장비를 조심스럽게 뒤로 기울이십시오.
 - b. 7/64 육각 키를 사용하여, 덮개를 장비 바닥판에 고정하고 있는 나사 4개를 푸십시오.



c. 장비를 피트에 조심스럽게 뒤로 기울이십시오.기기를 뒤로 젖히면 4개 나사가 빠집니다. 잃어버리지 않도록 주의하십시오.

d. 덮개를 바닥판에서 똑바로 들어 올리십시오.



- 4. 장비 덮개를 교환하십시오.
 - a. 장비 바닥판의 제자리에 덮개를 조심스럽게 놓으십시오.
 - b. 장비가 열 핀에 놓이도록 장비를 뒤로 기울이십시오.
 - c. 이전에 분리한 4개 나사를 삽입하십시오.모든 나사를 조이십시오.
 - d. 장비를 피트에 뒤로 기울이십시오.
- 5. 크리스탈 플레이트를 다시 장착하십시오.
 - a. 플레이트를 플랜지 위의 제자리에 놓으십시오.
 - b. 와셔를 나사에 끼우고 세 구멍에 나사를 넣으십시오.나사를 구멍에 똑바로 내려야 합니다.
 - c. 크리스탈 플레이트를 누른 상태에서 플레이트 양쪽에 있는 2개의 나사를 조인 후 장비 뒷면에 가장 가까운 세 번째 나사를 조이십시오. 손가락으로 모든 나사를 조이십시오.

나사가 자유롭게 회전하는 경우 나사를 위로 올린 후 다시 삽입하여 정렬을 수정하십시 오.

6. 분광기를 켜십시오.

- a. 전원 코드를 장비에 연결하고 전원을 켜십시오.
- b. 퍼지 라인을 장비에 다시 연결하고, 적용되는 경우 퍼지를 켜십시오.
- 7. 분광기의 성능을 확인하십시오.
 - a. 분광기 광학장치를 조절하십시오.
 - i. OMNIC Paradigm 소프트웨어의 대시보드에서 Acquire Data(데이터 획득) > Diagnostics(진단) > Align Spectrometer(분광기 조절)로 이동하십시오.
 - ii. Start(시작)를 선택하여 조절을 시작하십시오.조절이 정상적으로 완료되면 화면에 'Align complete(조절 완료)'메시지가 표시됩니다.
 - b. 레이저를 보정하십시오
 - i. Acquire Data(데이터 획득) > Diagnostics(진단) > Laser Calibration(레이저 보정)으로 이동하십시오.
 - ii. Start(시작)를 선택하여 보정을 시작하십시오.

레이저 보정이 완료될 때까지 몇 분 정도 걸릴 수 있습니다. 보정이 완료되면 디스플 레이에 'Measurement complete(측정 완료)' 메시지가 표시되고, 레이저 주파수의 변화 가 표시됩니다.

장비가 열평형에 도달할 때까지 최대 12시간 정도 걸립니다.최고의 결과를 얻으려면 처음 12시간 작동 후 조절과 보정을 다시 실행하십시오.

- c. Nicolet Summit OA PV Test(Nicolet Summit OA PV 테스트) 작업 플로우를 실행하십 시오.성능 확인 작업 플로우에서는 기준 물질을 사용하여 일련의 표준 테스트를 실행하고 장비 작동을 확인하고 데이터의 정확성을 확보합니다.
 - i. 대시보드로 돌아가십시오.
 - ii. Workflows(작업 플로우) 창으로 스크롤 이동하여 Nicolet Summit OA PV Test (Nicolet Summit OA - PV 테스트)를 선택하십시오.
 - iii. 작업 플로우를 마우스 오른쪽 버튼으로 클릭하고 Run(실행)을 선택하여 작업 플로우 를 시작하십시오.

건조제 교환

OMNIC Paradigm 소프트웨어를 사용하여 장비 습도를 모니터링하십시오.필요한 대로 또는 조 직의 정책에 따라 건조제를 교환하십시오.

건조제를 교환하려면 장비 덮개를 탈거하고 건조제 카트리지를 분리하고 건조제 캐니스터와 O링을 교환해야 합니다.

필요한 도구

시작하기 전에, 다음 도구를 구비해야 합니다.

- 7/64 육각 키
- 건조제 교환 킷
- 3/32 육각 키
- #2 십자 드라이버
- 장갑, 손가락 골무 또는 실험용 티슈(습도 표시기 취급용)

경고

폭발 위험 방지.



장비 덮개를 탈거하고 건조제실을 열기 전에,장비 전원을 끄고,전원 코드 를 분리하고,시스템에서 샘플을 모두 제거하십시오.

가연성 액체 또는 가스가 건조제실에 들어가면 폭발 위험이 발생합니다.이 러한 경우 즉시 당사에 연락하고,이 조건이 해소될 때까지 장비의 전원을 켜지 마십시오.

알림

건조제 덮개를 분리했을 때 장비에 아무런 물체도 떨어지지 않도록 하십시오.



◆ 건조제를 교환하려면

- 1. 분광기를 끄십시오.
 - a. 분광기를 끄고, 전원 버튼을 누르십시오.
 - b. 전원 코드를 분리하십시오.
 - c. 퍼지를 끄고, 해당하는 경우 장비에 연결된 퍼지 라인을 분리하십시오.

참고 건조제 카트리지(퍼지 커넥터 포함 또는 제외)를 사용하여 필요한 습도를 유지 할 수 있습니다.아래 그림은 퍼지되지 않은 건조제 카트리지 버전입니다.건조제 캐 니스터 교환 지침은 두 카트리지에 동일합니다.

2. "크리스탈 플레이트 탈거"

- 3. 장비 덮개를 탈거하지 마십시오.
 - a. 장비가 열 핀에 놓이도록 장비를 조심스럽게 뒤로 기울이십시오.
 - b. 7/64 육각 키를 사용하여,덮개를 장비 바닥판에 고정하고 있는 나사 4개를 푸십시오.



c. 장비를 피트에 조심스럽게 뒤로 기울이십시오.기기를 뒤로 젖히면 4개 나사가 빠집니다. 잃어버리지 않도록 주의하십시오.

d. 덮개를 바닥판에서 똑바로 들어 올리십시오.



- 4. 건조제 카트리지를 분리하십시오.
 - a. #2 십자 드라이버를 사용하여 건조제 카트리지에서 두 계류 나사를 풀고, 카트리지를 수 직으로 들어 올려서 장비에서 꺼내십시오.



b. 큰 O 링을 분리하여 버리십시오.

Thermo Scientific



참고 O 링을 건조제실에 떨어뜨리지 않도록 주의하십시오.

5. 건조제 캐니스터를 분리하고 습도 표시기를 교환하십시오.



a. 0.05인치 육각 키를 사용하여 포화된 두 건조제 캐니스터를 건조제 카트리지에 고정하고 있는 8개 나사를 푸십시오(약 2 바퀴). 캐니스터를 분리하십시오.



b. 건조제 카트리지를 위쪽이 아래를 향하도록 뒤집고 창에서 습도 표시기를 벗기십시오. 사용한 표시기를 버리십시오.

알림

새 습도 표시기를 취급할 때 항상 실험실용 장갑 또는 손가락 골무를 착용하거나 실험 용 티슈를 사용하십시오. 피부의 오일이나 습기는 표시기를 변색시킬 수 있습니다.

- 6. 새 건조제 캐니스터를 설치하십시오.
 - a. 새 건조제가 들어 있는 밀봉된 패킷(밀봉하여 건조제를 유지해야 함)을 개봉하고 새 습 도 표시기를 제거하십시오.
 - b. 건조제 카트리지를 위쪽이 아래를 향하게 뒤집고 표시기의 편평한 가장자리가 카트리 지의 가장자리와 일치하도록 창에 대고 파란색 표시기를 누르십시오. 카트리지를 올바 르게 다시 뒤집으면 창을 통해 문자가 보여야 합니다.

알림

카트리지를 다시 설치할 때 표시기가 창에 단단히 밀착되어 건조제실에서 느슨해지 지 않도록 하십시오.

- c. 캐니스터를 건조제 카트리지의 채널에 삽입하고 카트리지 채널과 수평이 될 때까지 8개 나사를 조이십시오(약 2 바퀴).
- 7. 건조제 카트리지를 삽입하십시오.
 - a. 패킷에서 새 O 링을 분리하고 장비 바닥판에 설치하십시오.

새 O 링을 제자리에 눌러서(굴리지 않음) 홈에 올바르게 장착되도록 하십시오.

b. 작업자가 장비 앞에 서서 습도 표시기의 글자를 읽을 수 있도록 건조제 카트리지를 분광 기에 삽입하십시오.

카트리지가 O 링 위에 올바르게 장착되었는지 확인한 후 #2 십자 드라이버를 사용하여 두 계류 나사를 조이십시오.

- 8. 장비덮개를 교환하십시오.
 - a. 장비 바닥판의 제자리에 덮개를 조심스럽게 놓으십시오.
 - b. 장비가 열 핀에 놓이도록 장비를 뒤로 기울이십시오.
 - c. 이전에 분리한 4개 나사를 삽입하십시오.모든 나사를 조이십시오.

- d. 장비를 피트에 뒤로 기울이십시오.
- e. "크리스탈 플레이트 교환".
- 9. 분광기를 켜십시오.
 - a. 전원 코드를 장비에 연결하고 전원을 켜십시오.
 - b. 퍼지 라인을 장비에 다시 연결하고,적용되는 경우 퍼지를 켜십시오.
- 10. 분광기의 성능을 확인하십시오.
 - a. 분광기 광학장치를 조절하십시오.
 - i. OMNIC Paradigm 소프트웨어의 대시보드에서 Acquire Data(데이터 획득) > Diagnostics(진단) > Align Spectrometer(분광기 조절)로 이동하십시오.
 - ii. Start(시작)를 선택하여 조절을 시작하십시오.조절이 정상적으로 완료되면 화면에 'Align complete(조절 완료)'메시지가 표시됩니다.
 - b. 레이저를 보정하십시오
 - i. Acquire Data(데이터 획득) > Diagnostics(진단) > Laser Calibration(레이저 보정)으로 이동하십시오.
 - ii. Start(시작)를 선택하여 보정을 시작하십시오.

레이저 보정이 완료될 때까지 몇 분 정도 걸릴 수 있습니다. 보정이 완료되면 디스플 레이에 'Measurement complete(측정 완료)' 메시지가 표시되고, 레이저 주파수의 변화 가 표시됩니다.

장비가 열평형에 도달할 때까지 최대 12시간 정도 걸립니다.최고의 결과를 얻으려면 처음 12시간 작동 후 조절과 보정을 다시 실행하십시오.

- c. Nicolet Summit OA PV Test(Nicolet Summit OA PV 테스트) 작업 플로우를 실행하십 시오.성능 확인 작업 플로우에서는 기준 물질을 사용하여 일련의 표준 테스트를 실행하고 장비 작동을 확인하고 데이터의 정확성을 확보합니다.
 - i. 대시보드로 돌아가십시오.
 - ii. Workflows(작업 플로우) 창으로 스크롤 이동하여 Nicolet Summit OA PV Test (Nicolet Summit OA - PV 테스트)를 선택하십시오.
 - iii. 작업 플로우를 마우스 오른쪽 버튼으로 클릭하고 Run(실행)을 선택하여 작업 플로우 를 시작하십시오.

소스 교환

IR 소스는 분광기 하단에서 쉽게 액세스할 수 있으며, 분광기 덮개를 분리하지 않고도 교환할 수 있습니다.

필요한 도구와 품목

- #1 십자 드라이버
- 소스 교환 킷

주의



소스를 교환하기 전에,장비를 끄십시오.당사가 제공하는 교환 부 품으로만 소스를 교환하십시오.

◆ 소스를 교환하려면

- 1. 분광기를 끄고 분리하십시오.
 - a. 장비를 끄고 전원 코드를 분리하십시오.
 - b. 장비에서 모든 케이블을 분리하십시오(이더넷 케이블, USB 기기 등).
 - c. 분광기가 퍼지된 경우 분광기에서 퍼지 라인을 분리하십시오. 자세한 내용은 <u>"퍼지 킷</u> <u>설치 및 유지보수"</u> 항목을 참조하십시오.



- 2. 소스를 분리하십시오.
 - a. #1 십자 드라이버를 사용하여 소스를 제자리에 고정하고 있는 3개 계류 나사를 완전히 푸십시오.



b. 베일을 분리하십시오. 한 손으로 장비를 고정하고 베일을 세게 당겨서 장비에서 소스를 바로 꺼내십시오.



알림

절대 맨손가락으로 소스 요소에 접촉하지 마십시오. 요소에 피부 오일이나 다른 퇴적물이 있으면 수명이 짧아집니다. 소스 요소를 취급할 때 항상 깨끗한 손가락 보호 골무, 장갑 또 는 깨끗한 실험용 티슈를 사용하십시오.

- 3. 새 소스를 삽입하십시오.
 - a. 3개 계류 나사가 직선인지 확인한 후 소스 장착부가 장비와 수평이 될 때까지 소스를 캐 비티에 천천히 삽입하십시오.
 - b. 나사를 조여서 소스를 고정하고 베일을 보관 위치에 잠그십시오.
- 4. 분광기를 다시 연결하고 전원을 켜십시오.
 - a. 주의하여 장비를 수직 위치로 되돌리고 분리한 모든 케이블을 다시 연결하십시오.
 - b. 건조제를 교환하십시오. 지침은 "건조제 교환" 항목을 참조하십시오.

소스를 분리할 때 내부 광학장치가 주위에 노출되기 때문에,소스를 교환한 후 항상 건조제 를 교환해야 합니다.

경고	
	감전 위험 방지.
	접지 연결을 올바르게 하고 감전 위험을 방지하려면 도관 접지에 연결된 콘센트를 사용하지 마십시오. 접지는 주 배전함의 접지에 연결되어 있는 부도체 배선이어야 합니다.

- c. 전원 코드를 장비에 연결하고 전원 버튼을 눌러서 분광기를 켜십시오.
- 5. 분광기의 성능을 확인하십시오.
 - a. 분광기 광학장치를 조절하십시오.
 - i. OMNIC Paradigm 소프트웨어의 대시보드에서 Acquire Data(데이터 획득) > Diagnostics(진단) > Align Spectrometer(분광기 조절)로 이동하십시오.
 - ii. Start(시작)를 선택하여 조절을 시작하십시오.조절이 정상적으로 완료되면 화면에 'Align complete(조절 완료)'메시지가 표시됩니다.
 - b. 레이저를 보정하십시오
 - i. Acquire Data(데이터 획득) > Diagnostics(진단) > Laser Calibration(레이저 보정)으로 이동하십시오.
 - ii. Start(시작)를 선택하여 보정을 시작하십시오.

레이저 보정이 완료될 때까지 몇 분 정도 걸릴 수 있습니다. 보정이 완료되면 디스플 레이에 'Measurement complete(측정 완료)' 메시지가 표시되고, 레이저 주파수의 변화 가 표시됩니다.

장비가 열평형에 도달할 때까지 최대 12시간 정도 걸립니다.최고의 결과를 얻으려면 처음 12시간 작동 후 조절과 보정을 다시 실행하십시오.

- c. Nicolet Summit OA PV Test(Nicolet Summit OA PV 테스트) 작업 플로우를 실행하십 시오.성능 확인 작업 플로우에서는 기준 물질을 사용하여 일련의 표준 테스트를 실행하고 장비 작동을 확인하고 데이터의 정확성을 확보합니다.
 - i. 대시보드로 돌아가십시오.
 - ii. Workflows(작업 플로우) 창으로 스크롤 이동하여 Nicolet Summit OA PV Test (Nicolet Summit OA - PV 테스트)를 선택하십시오.

 iii. 작업 플로우를 마우스 오른쪽 버튼으로 클릭하고 Run(실행)을 선택하여 작업 플로우 를 시작하십시오. [별도 공란]