Nicolet Summit LITE

FTIR 光谱仪



用户手册

269-345500 修订版 C 2021 年 4 月

© 2020 Thermo Fisher Scientific Inc. 保留所有权利。

Thermo Fisher Scientific Inc. 向购买产品的客户提供本文档,用作产品操作指南。本文档 受版权保护,未经 Thermo Fisher Scientific Inc.书面许可,严禁复制本文档或本文档中的 任何部分。

本文档的内容可能随时更改, 恕不另行通知。本文档中的所有技术信息仅供参考。本文档中的系统配置和规格将取代购买者先前获得的所有信息。

Thermo Fisher Scientific Inc. 不保证本文档的完整性和准确性,而且对于可能因使用本 文档(即使是在正确遵循本文档中的说明信息的情况下)而导致的任何错误、疏忽、损害 或损失, Thermo Fisher Scientific Inc. 概不负责。

本文档不属于 Thermo Fisher Scientific Inc. 和购买者之间的销售合同的一部分。本文档不 应以任何方式约束或修改任何"销售条款和条件",若两份文档信息发生冲突,则以"销售 条款与条件"中的信息为准。

仅供研究使用。此仪器或附件不是医疗器械,因此不适合用于预防、诊断、护理或治疗 疾病。



目录

| 欢迎 | . 1 |
|-------|-----|
| 使用的约定 | . 2 |
| 定购零件 | . 3 |
| 联系我们 | . 3 |

| Nicolet Summit LITE 光谱仪入门 | 5 |
|----------------------------|----|
| Summit 光谱仪到货前的准备工作 | 6 |
| 使用 Summit 光谱仪的个人安全事项 | 8 |
| 拆箱并开始使用光谱仪 | |
| 安装采样附件 | |
| 了解光谱仪灯带 | 23 |
| 使用 OMNIC Anywhere 应用程序查看数据 | |
| | |

| 选配和附件 | 27 |
|---------------------|----|
| 适用于 Summit 光谱仪的采样附件 | |
| 安装和维护吹扫套件 | |

| 维护 | |
|----------------|----|
| Summit 光谱仪维护计划 | |
| 光谱仪光学元器件准直 | 45 |
| 校准激光频率 | 46 |
| 清洁光谱仪 | 47 |
| 更换干燥剂 | |
| 更换光源 | |
| 更换样品仓窗片 | |

第1章

欢迎

Thermo Scientific Nicolet Summit LITE 傅里叶变换红外 (FTIR)光谱仪可通过各种附件采集中红外 光谱范围中的数据,针对样品材料执行化学分析。该系统集成了验证功能、强大的软件套件以 及其他众多功能,显著简化了数据采集。用户可自行安装选配硬件并执行多种维修和维护程 序。本文档或其他随附文档包含用户所需信息。

请务必首先阅读系统随附的安全指南,然后再操作仪器。

注意 建议仪器始终保持密封和干燥并/或对仪器进行吹扫。因未保持密封和干燥和/或未进 行吹扫而导致的设备损坏不在保修范围内。如对此要求有疑问,请联系我们。

使用的约定

安全防范措施和其他重要信息使用以下格式:



注意 包含有帮助的补充信息。

Thermo Scientific

定购零件

如需订购部件,请联系我们。

如果需要将仪器或附件送修,请先致电或发邮件联系我们,了解运送要求或其他说明。

联系我们

如需技术支持,请联系:www.thermofisher.com

[本页有意留为空白]

第2章

Nicolet Summit LITE 光谱仪入

为了顺利使用新的 Nicolet Summit LITE 光谱仪,请在仪器到货前阅读场地和安全要求。仪器到货后,需拆开包装、安装仪器并连接其他附件。

| Summit 光谱仪到货前的准备工作 | 6 |
|----------------------------|----|
| 使用 Summit 光谱仪的个人安全事项 | 8 |
| 拆箱并开始使用光谱仪 | |
| 安装采样附件 | |
| 了解光谱仪灯带 | 23 |
| 使用 OMNIC Anywhere 应用程序查看数据 | 26 |

Summit 光谱仪到货前的准备工作

在光谱仪到货前,需要为其准备一个合适的工作区,确保获得最精准的数据及持久耐用的最佳性能。

本文介绍光谱仪到货时需要执行的操作,以及可能影响光谱仪性能的环境和电气因素。如需更加全面地了解这些要求,请参阅光谱仪的场地和安全信息。

光谱仪拆箱

在光谱仪到货后,打开包装箱前需要立即执行两大重要步骤:

• 检查装运箱外部是否损坏

如果发现任何损坏迹象,请联系我们或当地分销商以获取相关指示。

•静置光谱仪,使其温度达到室温

在装运箱内,光谱仪密封在一个塑料袋中以保持干燥。**打开密封塑料袋前,必须静置 24 小时,等待光谱仪的温度达到室温。**如果在光谱仪温度达到室温前就将塑料袋打开,会形成 冷凝水,损坏内部光学组件,导致永久损坏。

保修范围不包括因移动方法不当或在仪器达到室温前拆除密封塑料袋所导致的损坏。

准备工作区

在仪器到货前,确保工作区能够正确安放光谱仪。除为光谱仪预留空间外,还需要考虑多项环 境和电气要求。

光谱仪尺寸

Summit 光谱仪需要的空间相当小,但仍应在仪器周围预留一定空间,以供通风孔散热并且便于操作光谱仪的端口、电源开关和电缆。

- Summit 仪器重量:9.6千克(21磅)
- 尺寸(宽 x高 x 深):34 厘米 x 24 厘米 x 32 厘米(13.3 英寸 x 9.6 英寸 x 12.7 英寸)

环境因素

Summit 光谱仪是一款坚固耐用的仪器,可在很多环境中使用,但其在相对无尘和低湿度环境中能够达到最佳性能。当温度介于 15 和 35 °C 之间时,光谱仪可稳定运行;在温度处于 20 到 22 °C 之间时,其性能达到最佳状态。

Thermo Scientific

潮湿会导致仪器内部形成冷凝水,损坏内部组件。通常可以采取几种预防措施防止仪器受潮:

- 保留仪器的干燥剂,在仪器存放期间也同样如此。
- 避免温度快速变化。
- 使仪器远离冷热空气源,例如靠近暖气和空调通风口或者大窗户。
 如果光谱仪存放在特别潮湿的环境中,应考虑安装吹扫气体套件。

电气要求

光谱仪必须使用专用的不间断电源供电,不能存在以下问题:

- 电压中断
- 瞬态尖峰
- 频移
- 其他线路干扰

如果怀疑电源有问题,建议对电源质量进行审核。有关详细信息,请联系我们或您所在地的电力局。

电气服务规格

下表列出了电气服务规格。如果您对这些要求有任何问题,请联系我们在当地的服务代表。

| 要求 | 规格 |
|------|---------------------------------|
| 输入电流 | 最大 1.5 A |
| 输入电压 | 100 至 240 VAC |
| 线路频率 | 50至60次 |
| 线路干扰 | 电压突降、电涌或其他线路干扰不得超出输入电压的 10% |
| 噪声 | 小于 2 V(普通模式) 小于 20 V(正常模式) |

Thermo Scientific

使用 Summit 光谱仪的个人安全事项

虽然 Nicolet Summit LITE 光谱仪采用保障安全性的仪器设计方案,但用户仍需采取一些措施预防正常使用和维护过程中潜在的危险。

注意事项



本手册将介绍用户需要了解的潜在危险,但并非面面俱到。使用仪器之前,请参阅光谱仪的场地和安全信息,查看有关这些潜在危险的完整说明。

正常使用期间的潜在危险

正常使用时的主要危险来源包括:

- 潜在危险样品和溶剂
- 接触硒化锌 (ZnSe)
- 光谱仪中温度较高或散发热量的部件
- 仪器的激光器

了解潜在的危险并采取适当的预防措施,有助于避免人身伤害或仪器受损。

危害性样品和溶剂

如果正在使用或计划测量潜在危害性的样品或溶剂(例如压缩气体或腐蚀性/易燃性溶剂),请 采取专门的预防措施。

适当通风

光谱仪对通风无特殊要求,但对于特定类型的分析,可能需要额外进行通风。如需分析剧毒样品、将样品溶解在可能与红外光源相互作用的溶剂中或采集易燃气体时,请确保通风良好。

含卤代烃的溶剂的热解可产生盐酸(HCI)、氢氟酸(HF)或光气(COCl₂。

警告



避免吸入毒物。氢氯酸、氢氟酸和光气都是高毒性的。如果使用含 有卤化烃的溶剂,请确保工作区域通风良好。

挥发性和易燃性溶剂

光谱仪内部的红外线光源可能点燃具备易燃性和挥发性的样品和溶剂。处理易燃样品和溶剂 时,请采取以下措施:

- 在安装样品室窗片的情况下工作。
- 使用不产生火花和其他火源的有效通风系统确保工作区通风良好,预防可燃蒸气在仪器周围的空气中聚集。
- 请勿将可燃性溶剂或样品置于仪器附近。
- 不需要时,请勿将可燃性溶剂或样品长时间放在样品室。
- 使用清洁、干燥的空气或氮气吹扫光谱仪。

腐蚀性溶剂

在样品室内使用产生 HCI或 HF 的样品可能会严重损害系统。如果使用含有卤化物的溶剂,请使用清洁、干燥的空气或氮气吹扫仪器。

注意

因不吹扫仪器而造成的设备损坏不包括在保修范围内。

HCI和 HF 气体也可能损害样品室 KBr 窗片的涂层。如果计划经常使用腐蚀性溶剂,样品室请考虑换用 ZnSe 窗片。

生物危害或放射性物质和传染性物质

生物样品,例如组织、体液、传染性物质和血液),均存在传播传染病的风险。使用存在传染性风险的物质时,请务必遵循贵机构的"生物安全计划"方案。

有毒物质

Nicolet Summit OA 光谱仪包括硒化锌 (ZnSe) ATR 晶体, 其他 Summit 仪器可使用选配的 ZnSe 样品室窗口。





避免吸入和摄入毒物。

硒化锌 (ZnSe) 有毒。有关处理和接触控制的信息,请访问 www.specac.com 参阅制造商的 ZnSe 安全数据表。

热源

在正常使用期间,光谱仪的某些部件可能变得烫手。请注意光谱仪的红外线光源和通风孔附近。

图2-1:注意光谱仪的红外线光源和通风孔附近



红外光源的外表面位于仪器底部,可以变得非常烫手。请勿在操作过程中或刚使用完仪器时触 摸仪器底部。

光谱仪的通风孔位于仪器左侧。在正常使用期间,光谱仪通过这些通风孔将内部组件产生的热气散发到光谱仪外部。请在仪器通风孔附近预留足够的空间以便散热。

激光和光学安全

在正常使用光谱仪时,用户永远不会遭受超出安全等级的激光辐射。如果在维修过程中取下盖板,可能需要采取特殊的预防措施,例如使用护目镜。维修人员将在必要时向用户发送通知。

警告



避免人身伤害。

切勿凝视激光束或其反射光。即使更换发生故障的激光器,也不要随意触碰激光器,以免受到激光或高压的伤害。

维护期间的潜在危险

维护仪器时的潜在危险与正常使用期间的危险不同。在维护期间,主要的危险来源于吹扫仪器 和操作仪器内部组件。

吹扫仪器。

在湿度偏高的环境中,我们建议安装清洁、干燥的空气或氮气源,用于吹扫光谱仪。吹扫仪器 有助于保护其内部光学组件免受潮湿环境或腐蚀性溶剂的损伤,确保获得精度更高的结果。

有关购买和安装光谱仪吹扫套件的详细信息,请参阅"安装和维护吹扫套件"。



操作内部组件

通常来说,用户不应取下仪器盖或试图更换内部组件。然而,如需取下盖板进行维护,请务必 注意电击、灼伤和激光辐射的危险。



注意事项



避免灼伤危险。

内部组件在正常运行时可能变得非常烫手,尤其是红外线光源。更换组件 之前务必关闭仪器并至少等待 10 分钟。

波在人自人



避免人身伤害。

- 切勿凝视激光束或其反射光。切勿随意触碰激光器,以免受到激光或高压的伤害。
- •如果调节激光器或未按用户手册所述的步骤进行操作,用户可能遭受有害辐射。

更换干燥剂

警告

打开干燥剂室时,必须避免易燃液体或气体进入干燥剂室。有关更换干燥剂的说明,请参阅<u>"更</u>换干燥剂"。

危险



避免爆炸的发生。

在打开干燥剂室之前,请关闭仪器电源,拔下电源线,并取下系统上所有的 附件和样品。如果易燃液体或气体进入干燥剂室,可能导致爆炸。如果发生 以上情况,请立即联系我们,并且在事故处理之前不要接通仪器电源。

更换红外线光源

红外线光源在使用过程中变得非常烫手。如需更换红外线光源,请注意灼伤和爆炸危险。

| 注意事项 | |
|------|--|
| | 避免灼伤危险。 |
| | 光源在使用过程中变得非常烫手。关闭光谱仪后,务必等待其冷却 10分钟 以上,才能操作光源。 |

Thermo Scientific

危险



避免爆炸的发生。

从光谱仪取下光源前,请关闭仪器电源,拔下电源线,断开所有吹扫管线并 取下系统上所有的附件和样品。如果易燃液体或气体进入光源室,可能导 致爆炸。如果发生以上情况,请立即联系我们,并且在事故处理之前不要接 通仪器电源。

总结

Nicolet Summit LITE 光谱仪是一款安全耐用的仪器,但在使用和维护过程中存在一定危险性。 在正常操作期间,处理潜在危险性的样品和溶剂时需要格外小心,同时避免接触温度较高或散 发热量的光谱仪部件。在维护过程中,请采取适当的预防措施,避免在吹扫仪器、操作内部组 件或更换干燥剂时造成人身伤害或设备损坏。

拆箱并开始使用光谱仪

拆箱并设置 Thermo Scientific Nicolet Summit LITE 光谱仪时,需要将仪器放置在工作区,连接光谱仪,接通电源并进行简短的性能测试。

准备工作

注意

请等待24小时再将光谱仪从塑料包装袋中取出,避免形成冷凝水而损坏内部光学元器件。 因在仪器达到室温前打开密封塑料包装袋而造成的损坏不在保修范围内。

提醒:查看光谱仪的现场和安全信息,确保工作区安全无虞,环境条件适合使用仪器。

开箱和设置

光谱仪的开箱和设置步骤稍有不同,具体取决于购买的选件。

1. 拆开装运箱

将光谱仪从装运箱和包装袋中取出,然后拆除其他包装材料。

注意

抬起光谱仪时请小心谨慎。

• 拆开仪器包装时,请勿触摸样品室窗片。触摸窗片可能造成永久性损坏。



2. (可选)更换样品室窗片。

如果购买了 ZnSe 窗片选件,可以更换窗片。有关更换窗片的说明,请参阅<u>"更换样品仓窗片</u>"。

3. 开启光谱仪。

- a. 将电源与光谱仪相连, 然后连接正确接地的交流电源。
- b. 按下仪器的电源开关打开光谱仪。仪器预热时,光谱仪的灯条循环发出绿光;当仪器准 备就绪时,灯条呈绿色常亮状态。



- 4. 安装 OMNIC Paradigm 软件。
 - a. 插入 OMNIC Paradigm 软件 DVD。
 - b. 如果 DVD 未自动启动,请打开 DVD 文件然后单击"开始"以开始安装。按照屏幕上的提示安装 OMNIC Paradigm 软件。

软件安装最多可能需要20分钟才能完成。您将需要重新启动计算机来完成安装。

- 5. 连接光谱仪
 - a. 使用 USB A 型或 B 型电缆将光谱仪连接到计算机。光谱仪配有 USB B 型插座。
 - b. 打开 OMNIC Paradigm 软件。首次打开该应用程序时, 它将在"工作站"模式下打开。

| Open | nd Sample | Settings C | eshboard S | MÊ Ê | T Advanced ATF | Search | Check Out | O Intify Add | f To Library | | |
|------------------------|-------------------------------|------------|--------------|-----------------|----------------|--------|------------|------------------------|--------------|------|--|
| Measurements | Past week | ~ | | | | Search | 0 Ø | | | | |
| Measurement Name | | Last | Modified | Туре | | | | | | | |
| Workflows 🗸 | | | | | | | freise | | | More | |
| Name | | Date | reated | Last Modified | | | PTENDA | | | | |
| ATR Accessory - PHEL | R Qualification | 8/10/5 | 020 12:00 PM | 8/10/2020 12:00 | | th | orr | nc | | | |
| Nicolet FTIR - CP Qua | ification | 8/10/0 | 020 12:00 PM | 8/10/2020 12:00 | | u | | IIC | / | | |
| Nicolet FTIR - Factory | Qualification | 8/10/2 | 020 12:00 PM | 8/10/2020 12:00 | | 50 | nair | tifi | \sim | | |
| Nicolet FTIR - JP Qual | fication | 8/10/2 | 020 12:00 PM | 8/10/2020 12:00 | | 00 | | un | | | |
| Nicolet FTIR - PHEUR | Qualification | 8/10/2 | 020 12:00 PM | 8/10/2020 12:00 | | | | | | | |

c. 连接光谱仪后,OMNIC Paradigm 软件会自动识别光谱仪并在屏幕的右上角显示仪器类型。如果仪器未自动连接,则可能需要手动连接。

要手动连接,请转到"配置>连接性",然后从"连接性"对话框中选择 Summit LITE。

软件可能需要几分钟才能启动并连接到仪器。软件将会在屏幕的右上角显示连接状态。 当状态显示您已连接时,请继续执行下一步。 **注意** 如果在连接仪器之前打开软件中的任何菜单,则"准直"和"校准"将被禁用。若要 启用它们,请在连接仪器后单击工具栏中的"仪表板"。

6. 验证光谱仪性能。

执行以下步骤之前,请确保采样室内无采样附件。

- a. 光谱仪光学元器件准直
 - i. 从 OMNIC Paradigm 软件的仪表板,转到采集数据 > 诊断 > 光谱仪准直。

ii. 选择**启动**开始进行准直。成功完成准直后,屏幕将显示"准直完成"。

b. 校准激光。

- i. 转到采集数据 > 诊断 > 激光校准。
- ii. 选择启动开始进行校准。

激光校准可能需要几分钟才能完成。校准完成后,显示屏将显示"测量完成"并指示激 光频率变化。

仪器最多需要 12个小时才能达到热平衡。为获得最佳结果,请在首 12小时的运行后重新运行准直和校准。

- c. 运行 Nicolet FTIR 出厂验证工作流。性能验证和鉴定工作流使用光谱仪内部的聚苯乙 烯参比物进行一系列标准测试,验证仪器的运行情况并确保数据精度。
 - i. 返回仪表板。
 - ii. 滚动至工作流窗格,然后选择 Nicolet FTIR 出厂验证。
 - iii. 双击该工作流并选择运行启动工作流。
- 7. 插入采样附件。

如需安装选配的 Everest ATR 附件或 iD1 透射附件,请握住附件的前后手柄,将其放到光谱 仪上。采样附件安装在光谱仪底板上的两个定位销上方,通过强磁铁固定到位。

Nicolet Summit LITE 光谱仪入门



• 有关安装附件的详细信息,请参阅"安装采样附件"。

后续步骤

恭喜您! 仪器准备就绪,可以使用。有关使用光谱仪识别和分析样品的更多信息,请参阅 thermofisher.com/ftir-help 所提供的 OMNIC Paradigm 指南和教程。

安装采样附件

Nicolet Summit LITE 光谱仪可安装多种采样附件,包括Thermo智能附件以及需要额外安装适配器的附件。

智能附件和底板易于安装和拆卸,并可通过 OMNIC Paradigm 软件自动识别。

安装智能附件

Thermo Scientific[™] Everest[™] ATR 附件等智能附件安装在光谱仪底板的两个定位销上方,通过磁铁固定到位。

图2-2: Nicolet Summit 上显示的定位销和磁铁的位置。



1.定位销

2.磁铁

3.螺纹孔

- ◆ 安装智能附件
- 1. 握住附件的前后手柄,将其放入光谱仪样品室。磁铁可将附件固定到位。 图2-3:在 Nicolet Summit PRO上,将附件降下到位。



2. 如需对仪器进行吹扫,请将标有"连接附件"的吹扫管线连接到附件背面的吹扫入口。详细信息,请参阅<u>"安装和维护吹扫套件"</u>。



注意如需拆除附件,可在必要时断开吹扫管线连接,然后握住附件手柄并将其提起。不使用附件时,将其存放于柜子或箱子等无尘环境中。

Nicolet Summit LITE 光谱仪入门

安装底板固定附件

iD Base 适配器适用于多种FTIR取样附件,包括大型透射附件和其他专用附件。



注意 如果不连接附件,请勿安装 iD Base 适配器。强磁铁将适配器固定到位;可能需要用工具将其撬下。

如果slide-mounted(滑动固定)附件的尺寸对于 iD1 透射附件而言过大,可使用固定 iD 底座的滑动固定装置。



iD Base 适配器配有两组螺丝孔,因此可朝任一方向安装滑动固定装置。如需安装滑动固定装置,请将其放置在螺丝孔上方,然后插入并拧紧两个垫圈和螺丝。

使用高度调节螺丝升高或降低样品或附件,使其与红外光束对齐。



- ◆ 使用 iD Base 适配器安装附件
- 1. 将附件与 iD Base 适配器相连。
- 2. 将 iD Base 适配器放置于光谱仪底板中的两个定位销上。强磁铁可将适配器和附件固定到 位。
- 3. 一些专用的底板固定附件使用附件随附的4枚螺丝固定在仪器上。

安装滑动固定附件

iD1透射附件设计用于容纳气体或液体透射样品池以及具有2"X3"滑动底座的薄膜或颗粒支架。



- ◆ 安装滑动固定附件
- 1. 将附件滑入 iD1 透射附件中的一对插槽。

选择将附件中心置于光束焦点处的插槽中(由附件外壳上的箭头指定)。



了解光谱仪灯带

Thermo Scientific Nicolet Summit LITE 光谱仪配备的灯带可快速直观地指示产品质量通过/失败 以及仪器状态,十分方便快捷。

下表介绍了所有灯带信号。

预热中



智能采集

| 信号 | 系统状态 | 说明 |
|---------|-------------------------|-------------|
| 蓝灯呈脉冲状态 | 正在采集智能 背景 | 所有灯带呈蓝灯闪烁状态 |
| 蓝灯零星亮起 | 正在采集样品 数据(由用户发 起) | 一小部分蓝灯反复亮起 |

分析结果

| 信号 | 系统状态 | 说明 |
|-------|----------------------------------|---|
| 绿灯占比 | 匹配值或 QCheck 结果高 于阈值 | 绿灯占比可反映匹配值。例如, 占比约 90% 代表匹配值为 90。 |
| 橙色灯占比 | 匹配值或 QCheck 结果低 于阈值 | 橙色灯占比可反映匹配值。例 如,占比约 30%代表匹配值为 30。 |

诊断

| 信号 | 系统状态 | 说明 |
|------|------|---|
| 红灯闪烁 | 系统错误 | 整个灯带呈红色闪烁状态。有关 错误说明,请参阅 OMNIC Paradigm 软件中的 系统状态 。 |

使用 OMNIC Anywhere 应用程序查看数据

OMNIC Anywhere 是基于云技术的应用程序,可用于查看、浏览或共享所连接的电脑、苹果电脑、Andorid 或 iOS 设备中的数据。

图 2-4: Web 浏览器中的 OMNIC Anywhere



借助 Nicolet Summit LITE 光谱仪和免费 Connect 帐户,可在教室或实验室测量样品、将数据上 传到 Connect 帐户,并在宿舍或工作区的其他设备上查看、浏览或共享数据。

有关通过 OMNIC Anywhere 创建 Connect 帐户和查看数据的说明,请访问"FTIR 云课堂"。

第3章

选配和附件

Nicolet Summit LITE 可与选配的吹扫套件配合使用,同时兼容多种采样附件。

| 适用于 Summit 光谱仪的采样附件 | 28 |
|---------------------|----|
| 安装和维护吹扫套件 | 31 |

适用于 Summit 光谱仪的采样附件

Thermo Scientific Nicolet Summit LITE FTIR 光谱仪兼容 Thermo Scientific[™] Nicolet[™] iS5 iD 采样附件及数百种第三方开发的附件。

有关安装采样附件的说明,请参阅"适用于 Summit 光谱仪的采样附件"。

借助以下附件和扩展底座, Summit光谱仪可快速便捷地测量多种样品。

EVEREST ATR 附件



Everest[™] ATR 附件属于高性能、通用型的单次衰减全反射附件,非常适合分析液体、固体、糊状或粉末物质。Everest ATR 附件专为 Summit 光谱仪而设计。

<u>联系我们</u>,了解详细信息。

ID1透射附件



iD1透射附件便于对薄膜、颗粒、液体和气体进行采样,适用于多种透射附件。样品仓最大可装入 10 厘米的气体样品池和比色皿。

<u>联系我们</u>,了解详细信息。

Thermo Scientific

选配和附件

ID5/ID7 ATR 附件



iD5 和 iD7 ATR 附件专为 Nicolet iS[™]5 光谱仪设计,属于 Everest ATR 附件的替代产品。此类 ATR 附件为性能卓越的多用途附件,而 Everest ATR 附件则建议应用于 Summit 光谱仪。

有关详细信息,请参阅<u>"适用于 Nicolet[™] iS5 光谱仪的 iD5 ATR 附件"</u>或<u>"适用于 Nicolet[™] iS5 光谱</u> 仪的 iD7 ATR 附件"。

IDBASE扩展底座



iD 扩展底座兼容多种 第三方FTIR 采样附件,包括大型透射附件和其他专用附件。将 iD 拓展底座与样品支架配合使用,以获得更多选择。

详细信息,请参阅"iD Base 适配器"。

样品支架



将样品支架与 iD拓展底座配合使用,以适配尺寸对于 iD1 透射附件过大的样品支架附件。 联系我们,了解详细信息。

GOLDEN GATE ATR 附件



Golden Gate ATR 附件采用的蓝宝石砧可提供高压接点,同时配有坚固耐用的外壳,可用于分析从单一颗粒、纤维、坚硬固体到腐蚀性液体的多种样品。Golden Gate ATR 附件是进行常量采样的理想之选。

详细信息,请参阅"适用于 Nicolet iS5 FTIR 光谱仪的 Thermo Scientific™ Golden Gate ATR"。

安装和维护吹扫套件

使用干燥的空气或氮气吹扫光谱仪可防止内部组件受湿气和其他环境污染物的破坏。安装吹扫套件时,需要组装阀门和调节器、更换干燥剂盒、连接吹扫气体并设置压力和流速。

警告 避免爆炸的发生 切勿使用易燃、可燃或有毒气体吹扫此仪器。吹扫气不得含有油及其他活性物质。来自光源或激光吸收的热量可能会点燃吹扫气体中的易燃气体或活性物质。仅使用干燥的空气或氮气吹扫您的仪器。

注意

建议仪器始终保持密封和干燥并/或对仪器进行吹扫。因未正确干燥或吹扫仪器而造成的损坏不在保修范围内。如对此要求有疑问,请联系我们。

所需设置

安装吹扫套件之前,需要一个符合 Thermo Scientific Nicolet™Summit 光谱仪规范的干燥空气或 氮气源。有关详细信息,请参阅仪器的"场地和安全指南"。

注意为了达到最佳效果,吹扫气体应干燥到-70°C (-94°F)或更低的露点温度。

所需材料

吹扫套件包括:

- 双区域吹扫歧管
- 吹扫干燥剂组件,带有吹扫干燥剂盒和 O型环

除吹扫套件外,还需要以下工具和材料:

- 3/4 英寸开口扳手
- 11/16 英寸开口扳手
- •2号十字头螺丝刀
- 螺纹密封胶带("生胶带"或"特氟龙胶带")

◆ 安装吹扫套件

- 1. 将管道组件安装到吹扫气源。
 - a. 在吹扫气源上安装调节阀以及 1/4 英寸外螺纹管道接头或 3/8 英寸内螺纹管道接头。(选择适合吹扫气源的阀门和管道接头。)



b. 如果使用 1/4 英寸外螺纹管道接头,请继续执行下一步。

如果在吹扫气源上使用 3/8 英寸内螺纹管道接头,请安装随吹扫套件附带的 3/8 英寸转 1/4 英寸异径管接头。安装异径管接头前,使用接头密封胶带将其缠好,然后使用 11/16 英寸开口扳手紧固连接处。



c. 用接头密封胶带缠好异径管接头或 1/4 英寸外螺纹管道接头, 然后安装压力联轴器。使用 3/4 英寸开口扳手紧固连接处。



d. 将墙壁管道组件的外螺纹入口牢固地卡入快卸接头。



- 1.连接附件
- 2.连接光谱仪
- 3.流量计
- 4.调压器

5.快速接头

6.墙壁管道组件

- 2. 更换干燥剂盒。
 - a. 请关闭仪器电源,拔下电源线并取下系统上所有的附件和样品。



b. 拧下标准干燥剂盒上的两枚十字螺丝,然后向上提起干燥剂盒,将其从仪器中取出。



注意 如果标准干燥剂盒存放在干燥盒子或密封袋中(可使用吹扫干燥剂盒的包装袋),则可重复使用该干燥剂盒。

c. 拆掉大的 O型环并将其丢弃。



- d. 打开吹扫套件和密封干燥剂包装。
- e. 从包装中取出新的O型环,将其安装在仪器的基板上(参阅上图)。将新的O型环按压 (不要卷动)到位,确保正确安放到凹槽中。
- f. 从包装中取出干燥剂盒及吹扫管道接头,将干燥剂盒插入干燥剂舱中,确保其完全压在 O型环上,然后用两枚螺丝将其固定。

请注意干燥剂盒的方向是否正确。干燥剂盒安装完毕后,应当可从仪器正面看到干燥剂 盒上的标签。



1.吹扫入口 2.释压阀

3.吹扫干燥剂盒

- 3. 将吹扫气源连接到仪器。
 - a. 将墙壁管道组件的气流联轴器(标有"连接光谱仪")连接到干燥剂盒上的吹扫入口。
 - b. 安装之前卸下的所有采样附件。
 - c. 如需吹扫附件,请将管道组件的直流联轴器(标有"连接附件")连接到所安装附件上的吹 扫入口。
- 4. 设置吹扫气体控制装置。
 - a. 打开主调节阀并按下文所述设置吹扫气体控制装置(有关详细信息,请参阅设置吹扫气 体控制装置)。如果不对附件进行吹扫,请将附件的调压器设为零。

| 硬件 | 压力 (psig) | 流速 (scfh) |
|---------------|------------------|-----------|
| Summit 光谱仪 | 5 | 1 |
| Everest 附件 | 5 | 4-10 |
| iD1 透射附件 | 5 | 4-10 |
| iD3 ATR 附件 | 5 | 4-10 |
| iD5 ATR 附件 | 5 | 4-10 |
| iD7 ATR 附件 | 5 | 4-10 |
| iD Foundation | 5 | 4-10 |

b. 将电源线插入仪器, 然后接通电源。

c. 等待 30-60 分钟, 使仪器充分进行吹扫。

设置吹扫气体控制装置

正确设置吹扫气体控制装置将防止光谱仪受潮,并且不会引起振动。 为了达到最佳吹扫效果,吹扫气体应干燥到 70°C (94°F)或更低的露点温度。

◆ 设置吹扫气体控制装置

- 1. 打开主阀门, 使吹扫气体开始流过调节器。
- 2. 调节光谱仪的调压器, 直至仪表指示压力为 5 psig (34 kPa)。



- 1.设置附件流量
- 2.设置附件压力
- 3.设置光谱仪流量
- 4.设置光谱仪压力
- 3. 将光谱仪的流量计设为1 scfh (0.47 l/min)。
- 4. 如果正在使用经过吹扫的附件,请按下文所述设置附件的吹扫气体控制装置。
 - 附件压力, 5 psig (34 kPa)
 - 附件流量: 4-10 scfh (1.9-4.7 l/min)

注意

流速超过建议值会引起振动,从而影响数据质量。

检查和更换吹扫气体过滤器

当吹扫过滤器变为黄色或褪色、或者受碎屑或外来颗粒污染,请更换。

注意

建议仪器始终保持密封和干燥并/或对仪器进行吹扫。因未保持密封和干燥和/或未进行吹扫 而导致的设备损坏不在保修范围内。如对此要求有疑问,请联系我们。

图3-1:吹扫滤网位于压力阀下面的塑料碗内。



注意 如需订购部件,请联系技术支持。

◆ 更换吹扫气体过滤器

1. 通过主阀门关闭吹扫气体。请勿调小流量计或调压器。

2. 取下盛放过滤器的塑料碗,然后取出过滤器。(塑料碗和过滤器均可用手拧下。)



- 3. 安装新过滤器,然后重新安装塑料碗。
- 4. 打开主阀门,开始向仪器输送吹扫气流,并验证吹扫气体流速是否正确。

[本页有意留为空白]

第4章

维护

Nicolet Summit LITE FTIR 光谱仪几乎无需保养,也不需要进行日常维护,但应进行基本养护,确保仪器持续高性能运行。

| Summit 光谱仪维护计划 | 42 |
|----------------|------|
| 光谱仪光学元器件准直 | . 45 |
| 校准激光频率 | . 46 |
| 清洁光谱仪 | . 47 |
| 更换干燥剂 | . 48 |
| 更换光源 | 54 |
| 更换样品仓窗片 | 58 |

Summit 光谱仪维护计划

Thermo Scientific Nicolet Summit LITE FTIR 光谱仪几乎无需保养,也不需要进行日常维护,但应进行基本养护,确保仪器持续高性能运行。请遵循本文中的指导原则。为了获得最佳性能,光谱仪可保持通电状态。

注意

静电会永久损坏光谱仪中的重要组件。为防止此类损坏,请遵循以下建议:

- 断电之前,触碰光谱仪的金属底座,释放可能积聚的所有静电。
- 在仪器中安装更换件之前,先将更换件存放在保护包装中。

每周维护

验证光谱仪性能

该OMNIC Paradigm 软件包含用于 Summit 光谱仪的认证和性能验证 (PV) 工作流。认证工作流属于用户最为认可的标准行业认证测试(例如欧洲药典、日本药典等)。PV 工作流运行一系列标准测试,验证仪器运行状况,确保数据精确。仪器提供了所需的全部参考标准,由OMNIC Paradigm 软件控制。

我们建议,至少每周运行一次 PV 工作流或您需要的认证工作流。

清洁光谱仪

严格根据建议清洁光谱仪。请参阅"清洁光谱仪"。

每月维护

检查湿度试纸

如果空气过于潮湿,很容易使光谱仪的光学组件受损。光谱仪密封保存,其组件通过两个吸湿的干燥剂进行保护。在 OMNIC Paradigm 软件中监控仪器湿度。

如果空气过于潮湿,很容易使光谱仪的光学组件受损。光谱仪密封保存,其组件通过两个吸湿的干燥剂进行保护。其中的湿度试纸用于检测光谱仪内部的湿度水平。

您也可以通过 OMNIC Paradigm 软件监控湿度。在桌面版界面中,单击仪器图标查看系统状态,或在触屏版界面中,选择"诊断"图标。"环境"部分将列出湿度水平。

图4-1:湿度试纸的位置



每月至少检查一次湿度试纸,需要时更换干燥剂(参阅下表),或者购买并安装吹扫套件。详细 信息,请参阅<u>"安装和维护吹扫套件"</u>。

| 湿度试纸状态 | 含义 | 措施 |
|--------|----------------------------------|----------------|
| 蓝色 | 干燥剂完全填充 | 无 |
| 浅蓝 | 干燥剂逐渐达到水分饱和 状态,无法再提供足够的 保护 | 更换干燥剂 |
| 粉色或白色 | 干燥剂过期 | 更换干燥剂和湿度 试纸 |

表1-湿度试纸状态和建议

详细信息,请参阅"更换干燥剂"。

注意

我们建议始终将光谱仪密封并保持干燥,或者使用干燥的空气或氮气吹扫。因未保持密封和 干燥和/或未进行吹扫而导致的设备损坏不在保修范围内。如对此要求有疑问,请联系我们。

检查吹扫气体滤网

如果使用氮气或者干燥空气对光谱仪进行吹扫,至少应每月检查一次吹扫滤网。

如果滤网变黄、变色或沾上了污染物或外界颗粒物,应进行更换。详细信息,请参阅<u>"检查和更</u>换吹扫气体过滤器"。

图4-2: 吹扫滤网位于压力阀下面的塑料碗内。



光谱仪光学元器件准直

如果验证工作流(例如 Nicolet FTIR - 出厂验证工作流)指示测试失败,请对光谱仪的光学元器件进行准直。对光谱仪准直可优化到达检测器的能量,最大限度增强检测器信号。如果性能或验证测试失败,请对光谱仪的光学元器件进行准直、校准激光频率并再次运行测试。如果测试仍然失败,请联系您当地的技术支持代表获得帮助。

◆ 对光谱仪的光学元器件进行准直(触屏版界面)

- 1. 在主屏幕上,选择"诊断"图标 [1]打开诊断视图。
- 2. 打开"准直"选项卡并选择准直。

系统自动对光谱仪进行准直,准直完成后出现提示。

- ◆ 对光谱仪的光学元器件进行准直(操作界面)
- 1. 浏览至管理员区,选择"诊断"图标 [200]打开诊断视图。
- 2. 打开"准直"选项卡并选择准直。

系统自动对光谱仪进行准直,准直完成后出现提示。

- ♦ 对光谱仪的光学元器件进行准直(桌面版界面)
- 1. 选择采集数据 > 诊断 > 光谱仪准直。
- 2. 在"光谱仪准直"对话框中,单击开始。

系统自动对光谱仪进行准直,准直完成后出现提示。

校准激光频率

如果验证工作流(例如 Nicolet FTIR - 出厂验证工作流)指示测试失败,请校准激光频率。如果测试失败,请先执行光谱仪准直,然后校准激光频率并再次运行测试。如果测试仍然失败,请联系您当地的技术支持代表获得帮助。

◆ 校准激光(触屏版界面)

- 1. 在主屏幕上,选择"诊断"图标 [1] 打开诊断视图。
- 2. 打开"激光校正"选项卡, 然后选择校准开始校准。

完成激光校准后,将出现一条消息,指示频率是否已更改并显示新的频率。

◆ 校准激光(操作界面)

- 1. 导航至管理员区,然后选择"诊断"图标 [1]打开诊断视图。
- 打开"激光校正"选项卡,然后选择校准开始校准。
 完成激光校准后,将出现一条消息,指示频率是否已更改并显示新的频率。
- ◆ 校准激光(桌面版界面)
- 1. 选择采集数据 > 诊断 > 激光校正。
- 2. 在"激光校正"对话框中,单击**开始**校准激光。

系统完成激光校准后,将出现一条消息,指示频率是否已更改并显示新的频率。

清洁光谱仪

在清洁 Thermo Scientific Nicolet Summit LITE 光谱仪前,请将其关闭并断开电源。

清洁 Summit 光谱仪

使用蘸有少量中性肥皂液的柔软洁净的抹布轻轻清洁仪器外部。

灰尘会积聚在仪器背面的电子组件外壳上,可能影响散热,进而影响电子组件的使用寿命。

如需清理仪器背面的灰尘,请使用压缩空气吹走灰尘。清理光谱仪背面的灰尘时,请勿使用任何液体。



注意

切勿使用强力清洁剂、溶剂、化学物质或研磨剂;这些都会对涂层造成损坏。请勿使液体接触 样品室中的任何窗口。

注意

窗片很容易刮划和损坏。请勿触摸窗片或尝试对其进行清洁。灰尘不会影响信号,但指纹会 导致仪器性能下降并会对镜面或窗片造成永久性损坏。如需清理镜面或窗片上的灰尘,请用 清洁、干燥的空气或氮气流将灰尘吹走。切勿让任何液体接触仪器中的窗口或光学元件。

更换干燥剂

如果仪器未配备吹扫套件,或者吹扫装置关闭,必须监测光谱仪内部的湿度水平。如果湿度试 纸变成粉红色(可能是浅粉红色或接近白色),请更换干燥剂。每次更换干燥剂时,都要同时更 换湿度试纸。

更换干燥剂时,需拆下干燥剂盒并更换干燥剂和O型环。

需要的工具

开始之前,需要准备以下工具:

- 干燥剂更换套件
- •一把 0.05 英寸的六角扳手
- 一把2号十字头螺丝刀
- 手套、指套或实验室用纸(处理湿度试纸)

警告



避免爆炸的发生。

打开干燥剂室前,请关闭仪器电源,拔下电源线,并取下系统上所有的附件 和样品。让易燃性液体或气体进入干燥剂室具有爆炸危险。如果发生以上 情况,请立即联系我们,并且在事故处理之前不要接通仪器电源。

注意

拆除干燥剂盒盖时,确保仪器内不落入任何东西。

警告



只能使用我们提供的部件更换干燥剂。

◆ 更换干燥剂

- 1. 关闭光谱仪。
 - a. 如需关闭光谱仪,请按下电源按钮。
 - b. 拔下电源线。
 - c. 关闭吹扫装置(如果适用)并断开吹扫管线与仪器和所有附件之间的连接。

注意可以使用任一干燥剂盒(带有或不带吹扫接头)保持所需的湿度。下面的几张图 中显示了不带吹扫装置的干燥剂盒。两个干燥剂盒更换干燥剂的指示相同。

- 2. 拆下干燥剂盒。
 - a. 拆下仪器上安装的所有附件,取出所有样品。
 - b. 使用2号十字头螺丝刀松开干燥剂盒上的两枚栓系螺丝,向上抬起干燥剂盒,将其从仪器中取出。



c. 拆掉大的 O型环并将其丢弃。



注意 小心不要让 O型环掉到干燥剂盒中。

3. 拆除干燥剂,更换湿度指示。



如果摄入干燥剂罐(水合金属硅铝酸盐分子筛)中的内容物,会危害健康。如果您要丢弃饱和的干燥剂罐,请确保妥善处理。

a. 使用 0.05 英寸的六角扳手, 拧松(大约拧两圈)将两个饱和的干燥剂固定在干燥剂盒上的八枚螺丝。拆下干燥剂。



1.干燥剂盒

2.拆下这些螺丝(每面四枚)

3.干燥剂

b. 将干燥剂盒倒过来,将旧的湿度试纸从窗上剥离下来。丢掉用过的试纸。

注意

处理新的湿度试纸时,始终要佩戴实验室手套或指套,或者使用实验室用纸。皮肤上的油或水分会使试纸变色。

- 4. 安装新的干燥剂。
 - a. 打开装有新干燥剂的密封包(必须密封保存干燥剂),拆下新的湿度试纸。
 - b. 将干燥剂倒过来,然后将蓝色试纸压到窗上,使试纸的平直边缘与盒子的边缘对齐。将 干燥剂盒的正面朝上时,应能通过窗口看到试纸上的文字。

注意

确保将试纸紧压到窗上,以便在重新安装干燥剂盒时,试纸不会在干燥剂仓内松动掉落。

- c. 将干燥剂插入干燥剂盒的通道中,拧紧八枚螺丝(约拧动2圈),直到与干燥剂盒管道平 齐。
- 5. 插入干燥剂盒。
 - a. 从包装中取出新的O型环,将其安装在仪器的基板上。将新的O型环按压(不要卷动)到位,确保正确安放到凹槽中。
 - b. 将干燥剂盒插入到光谱仪中,使站立在仪器前的人员能够看清湿度试纸上的文字。
 确保干燥剂盒正确安放在O型环上方,然后使用2号十字头螺丝刀拧紧两枚栓系螺丝。
- 6. 开启光谱仪。
 - a. 将电源线插入仪器, 然后接通电源。
 - b. 如果适用,将吹扫管线重新连接到仪器,然后打开吹扫装置。
 - c. 将光谱仪重新连接到计算机。
- 7. 验证光谱仪性能。

执行以下步骤之前,请确保采样室内无采样附件。

- a. 光谱仪光学元器件准直
 - i. 从 OMNIC Paradigm 软件的仪表板,转到采集数据 > 诊断 > 光谱仪准直。
 - ii. 选择启动开始进行准直。成功完成准直后,屏幕将显示"准直完成"。
- b. 校准激光。
 - i. 转到采集数据 > 诊断 > 激光校准。
 - ii. 选择启动开始进行校准。

激光校准可能需要几分钟才能完成。校准完成后,显示屏将显示"测量完成"并指示激 光频率变化。

仪器最多需要 12个小时才能达到热平衡。为获得最佳结果,请在首 12小时的运行后重新 运行准直和校准。

- c. 运行 Nicolet FTIR 出厂验证工作流。性能验证和鉴定工作流使用光谱仪内部的聚苯乙烯参比物进行一系列标准测试,验证仪器的运行情况并确保数据精度。
 - i. 返回仪表板。
 - ii. 滚动至工作流窗格,然后选择 Nicolet FTIR 出厂验证。
 - iii. 双击该工作流并选择运行启动工作流。

更换光源

从光谱仪底部可轻松更换红外光源,无需拆下光谱仪盖板即可进行更换。

所需工具和材料

- 一把 1 号十字头螺丝刀
- 光源更换套件

注意事项



更换光源之前,请将仪器关闭。只能使用我们提供的更换件更换光源。

◆ 更换光源

- 1. 关闭光谱仪并断开其连接。
 - a. 关闭仪器并拔下电源线。
 - b. 断开仪器上的任何电缆, 例如以太网电缆或 USB 设备。
 - c. 如果正在对光谱仪或安装的附件进行吹扫,则应将光谱仪或附件上的吹扫管线断开(快速连接配件自动阻断气流)。详细信息,请参阅<u>"安装和维护吹扫套件"</u>。
 - d. 从仪器上拆下安装的所有附件。

注意事项



避免灼伤危险。

在使用仪器时,红外线光源会变得非常烫手。关闭仪器后,至少等待 10分钟再执行下一步。

Thermo Scientific

- 2. 拆卸光源。
 - a. 小心将仪器后倾,直到停靠在电子组件外壳上。



b. 使用1号十字头螺丝刀将固定光源的三枚栓系螺丝完全松开。



c. 将拉手转出。一只手固定仪器,稳定地拉动拉手,从而将光源平直拉出仪器。



注意

切勿直接用手指触碰光源器件。如果器件沾上了皮肤油脂或其他沉积物,会缩短其使用 寿命。处理光源器件时,应始终佩戴干净的指套、手套,或者使用清洁的实验室用纸。

- 3. 插入新光源。
 - a. 确保三枚栓系螺丝方向平直, 然后缓慢将光源插入光源腔体中, 直到光源固定架与仪器 平齐。
 - b. 拧紧螺丝以固定光源,将拉手卡入原位。
- 4. 重新连接光谱仪并接通电源。
 - a. 小心地将仪器恢复到直立位置,重新连接拔掉的电缆。
 - b. 更换干燥剂。相关说明,请参阅"更换干燥剂"。

拆除光源时,内部光学元器件会暴露在空气中,因此在更换光源后,必须更换干燥剂。



c. 将电源线插入到仪器,按下电源按钮打开光谱仪。

注意 在对光谱仪进行准直并验证其性能之前,不要安装任何采样附件

5. 验证光谱仪性能。

执行以下步骤之前,请确保采样室内无采样附件。

- a. 光谱仪光学元器件准直
 - i. 从 OMNIC Paradigm 软件的仪表板,转到采集数据 > 诊断 > 光谱仪准直。
 - ii. 选择启动开始进行准直。成功完成准直后,屏幕将显示"准直完成"。
- b. 校准激光。
 - i. 转到采集数据 > 诊断 > 激光校准。
 - ii. 选择启动开始进行校准。

激光校准可能需要几分钟才能完成。校准完成后,显示屏将显示"测量完成"并指示激 光频率变化。

仪器最多需要 12个小时才能达到热平衡。为获得最佳结果,请在首 12小时的运行后重新运行准直和校准。

- - i. 返回仪表板。
 - ii. 滚动至工作流窗格,然后选择 Nicolet FTIR 出厂验证。
 - iii. 双击该工作流并选择运行启动工作流。
- 6. 更换先前拆下的所有附件。

Thermo Scientific Nicolet Summit LITE FTIR 光谱仪样品仓两侧均有窗片密封,可将仪器与水分和 其他污染物隔开,但允许红外光束进出样品仓。应始终安装窗片,即使用干燥气体或氮气吹扫 系统,也需要安装。



警告



避免火灾和爆炸的发生。

仪器内的红外线光源是一个起火源。如果正在使用挥发性溶剂,请准备通风橱或其他无火花和起火源的主动式通风系统,并防止易燃蒸汽在仪器周围空气环境中聚集。

光谱仪出厂时已安装溴化钾 (Kbr) 窗片,并在安装窗片的情况下运行了所有操作测试和性能测试。KBr 窗片具有吸湿性,但涂有防护涂层。在此光谱仪上,KBr 窗片提供了介于 8,000 与 350cm⁻¹ 之间的光谱范围,并且没有明显的光吸收线。

为了让光谱仪达到最佳性能,样品仓窗片必须清晰(不模糊)和洁净(没有灰尘和指印)。如果窗 片被污染,光谱仪可能无法通过性能和认证测试。可以从我们这里订购新的窗片,并按照下述 指示自行安装。 注意

请勿让液体接触窗片。 窗片很容易刮划和损坏。请勿触摸窗片或尝试对其进行清洁。灰尘不会影响信号,但指纹会导致仪器性能下降并会对镜面或窗片造成永久性损坏。如需清理镜面或窗片上的灰尘,请用清洁、干燥的空气或氮气流将灰尘吹走。(不要使用罐中的压缩空气,污染物可能导致器件损坏) 在准备将新的窗片插入光谱仪之前,应将其存放在保护包装中。 如果从光谱仪上拆下使用中的 KBr或 ZnSe 窗片,应立即放入提供的保护包装中(含干燥剂),然后将包装封牢。 KBr 窗片无色透明, ZnSe 窗片为黄色。 需要的工具: Nicolet Summit ZnSe(或 KBr) 窗片更换套件 套件中包含以下物品: ZnSe(或 KBr)更换窗片 (2)

- 窗片更换工具
- 用于存放不使用的窗片的保护包装(包括干燥剂)

• 只能使用我们提供的更换件更换样品仓窗片。

• 丁腈手套



◆ 更换样品仓窗片

1. 关闭光谱仪。

注意 如果仪器正在吹扫,可在更换窗片时保持吹扫装置打开,防止周围空气进入光谱 仪。

- 2. 拆下光谱仪上的所有采样附件。
- 3. 拆下最先安装的窗片。
 - a. 将窗片更换工具内缘上的三个凸条与最先安装的窗片的外缘槽口对齐。



注意 窗片很牢固。一只手稳定住仪器,另一只手拆卸窗片。

b. 牢牢按住工具,逆时针旋转以松开窗片。

继续松开窗片直至可以取下。将工具向上倾斜,即可取下窗片,因此不需要直接接触窗片。



注意

使用丁腈手套操作窗片,只能碰触边缘进行操作。(避免接触窗片表面,即使佩戴手套也应避免接触。)

- c. 如果窗片仍然可以使用,可拿着窗片的塑料环,将其小心放到提供的包装中(含干燥剂)。将包装存放在清洁干燥的环境中。
- 4. 安装新窗片。
 - a. 拿着新窗片的塑料外圈,小心放入窗片更换工具中,窗片螺纹朝上。
 - b. 旋转窗片, 直至其外缘中的三个凹口与工具内缘的凸条对齐。
 - c. 与拆卸步骤相反,将工具和窗片倾斜至恰好能将窗片放在光谱仪开口上方的角度。
 - d. 轻轻压住光谱仪侧壁,同时沿顺时针方向缓慢旋转工具,确保窗片螺纹正确啮合。

注意

开始时窗片应该容易转动。如果不是这样,请反向旋转,然后重新插入,以免破坏螺纹。

- e. 继续顺时针旋转工具,直至感觉到O型环压紧,然后再多转1/8圈。
- 5. 重复执行上述第3步到第4步,更换第二个窗片。
- 6. 打开仪器电源,连接至计算机,然后启动 OMNIC Paradigm 软件。
- 7. 验证光谱仪性能。

执行以下步骤之前,请确保采样室内无采样附件。

- a. 光谱仪光学元器件准直
 - i. 从 OMNIC Paradigm 软件的仪表板,转到采集数据 > 诊断 > 光谱仪准直。
 - ii. 选择启动开始进行准直。成功完成准直后,屏幕将显示"准直完成"。
- b. 校准激光。
 - i. 转到采集数据 > 诊断 > 激光校准。
 - ii. 选择启动开始进行校准。

激光校准可能需要几分钟才能完成。校准完成后,显示屏将显示"测量完成"并指示激 光频率变化。

仪器最多需要 12个小时才能达到热平衡。为获得最佳结果,请在首 12小时的运行后重新运行准直和校准。

- c. 运行 Nicolet FTIR 出厂验证工作流。性能验证和鉴定工作流使用光谱仪内部的聚苯乙 烯参比物进行一系列标准测试,验证仪器的运行情况并确保数据精度。
 - i. 返回仪表板。
 - ii. 滚动至工作流窗格,然后选择 Nicolet FTIR 出厂验证。
 - iii. 双击该工作流并选择运行启动工作流。
- 如果使用运行认证 (OQ) 测试跟踪仪器性能,建议在更换样品仓窗片后,重新运行所选 OQ 测试(特别是更改了窗片类型时)。有关详细信息,请参阅联机帮助中的"Summit 光谱仪认 证"一文。
- 9. 更换之前从光谱仪上拆除的附件。