

TLC Explorer 记录系统 1.52610.0001



用户操作手册



目录

1	简介.	•••••		6		
	1.1	TLC Ex	plorer的设计用途	6	2	
	1.2	安全与处理说明				
		1.2.1	在以下情况中,切勿操作TLC Explorer	7		
		1.2.2	安装要求	8		
		1.2.3	紫外线照射	8	2	
		1.2.4	一般操作注意事项	9	3	
		1.2.5	用户管理与信息安全	9		
	1.3	TLC简:	介	10		
		1.3.1	TLC工作原理	10	А	
		1.3.2	比移值R _f	11		
	1.4	关于本	手册	11		
2	了解T	LC Exp	lorer	12		
	2.1	TLC Ex	plorer的正面	12	5	
		2.1.1	· F1 - 电源按钮	12		
		2.1.2	F2, F3, F4 - 光照按钮和指示灯	13		
		2.1.3	F5, F6 - 抽屉和抽屉打开指示灯	13		
		2.1.4	F7, F8 - 检测窗和滑块	13	6	
	2.2	TLC Explorer的背面				
		2.2.1	B1 – 用于插接电源的USB-C端口	14		
		2.2.2	B2, B3 - USB-A端口	14		
		2.2.3	B4 - 以太网端口	14	7	
		2.2.4	B5 - WiFi天线连接器	15		
		2.2.5	B6 - 序号	15		
		2.2.6	B7 - 背面罩板	15		
		2.2.7	B8 - 光照装置	15	8	
	2.3	TLC Ex	plorer的底板	16		
		2.3.1	P1 - 底板区域	16		
		2.3.2	P2 - 边框	16		
		2.3.3	P3, P4 - 对准销和底座	16	9	
	2.4	附件…	·	17		
		2.4.1	A1 - WiFi天线	17		
		2.4.2	A2, A3 - 电源适配器和电源线	17		
		2.4.3		18	10	
3	初次何	刀次使用19				
	3.1	拆除包装与交付清单19				
	3.2	适合安全操作的设置				
	3.3	插电和	'首次查看'底板和色谱板	20	11	
	3.4	连接笔	记本电脑	21		
	3.5	创建首	位用户	23		
	3.6	关闭设	备	25		
					12	

4	操作	皆南	26
	4.1	典型工作流程	26
		4.1.1 主菜单	
	4.2	准备设备	
	4.3	目视检查	29
	4.4	捕捉图像	29
	4.5	加载现有图像集	
	4.6	选择和编辑色谱板	
	4.7	定义轨迹	35
	4.8	创建评估	37
	4.9	注解板图像	
	4.10	绘制和比较轨迹密度图	40
		4.10.1 绘制密度图	40
		4.10.2 添加斑点的名称或注释	41
	4.11	定量分析	43
		4.11.1 准备定量评估	43
		4.11.2 读取定量测量结果和微调	46
	4.12	导出图像和评估结果	47
		4.12.1 '导出原始裁切图像'(所有标签页)	48
		4.12.2 '导出显示的图像'(所有标签页)	48
		4.12.3 '导出注释'(仅 '注释')	48
		4.12.4 '导出斑点'('注释'和'比较')	49
		4.12.5 '导出比较'('比较'和 '定量')	49
		4.12.6 '导出定量'(仅 '定量')	49
		4.12.7 下载或保存导出的文件	49
	4.13	生成有关评估的报告	50
	4.14	管理文件和图像集	51
		4.14.1 导出图像集和下载图像	52
		4.14.2 复制或导入图像集	53
		4.14.3 删除图像集	53
	4.15	重置或关闭设备	53
5	用户管	管理和设备设置	54
	5.1	更改活动的用户	55
	5.2	添加和删除用户	55
	5.3	设备名和序号	56
	5.4	时间和日期	56
	5.5	数字光照校正	56
	5.6	杂散光检测	

目录

I		
•		
٠	•	

6	备份和	和数据迁移	
	6.1	生成备份	
	6.2	恢复备份	
	6.3	将用户数据迁移至不同的TLC Explorer58	
7	软件酮	更新	
	7.1	当前软件版本	
	7.2	获取软件更新	
	7.3	安装更新	4
	7.4	恢复前一软件版本	
	7.5	恢复出厂设置61	
8	连接1	۲LC Explorer	
	8.1	TLC Explorer自己的WiFi	
	8.2		
	8.3	集成到现有WiFi网络中64	
	8.4	重置网络连接64	
	8.5	将TLC Explorer作为网络位置(SMB)安装64	
9	故障排	非除65	
	9.1	识别错误65	
	9.2	可行的补救措施和解决方案66	
10) 清洁、	维护和停用	
	10.1	清洁	
	10.2	联系技术支持	
	10.3	订购附件和更换部件	
	10.4	更换光照装置	
	10.5	停用、储存和处置	
11	术语	表	
12	技术数	数据	
	12.1	设备测量和规格	
	12.2	光照频谱	

1.1 TLC Explorer的设计用途

TLC Explorer用于记录和评估薄层色谱板。它提供在完成TLC色谱板的实验室色谱展开工作流程后,运用目视检查和电子方法记录TLC色谱板图像的灵活、安全的快速方法。该设备包含三个不同的LED光源,它们在直接照射模式下从上方照射色谱板,适合检测不同的物质种类:

- · **白光**用于检测薄层色谱图上的可见有色物质。
- 长波紫外线(UV-A) (366 nm)一般用于激发荧光。在这种光照下,荧光物质和物体在黑暗的背景中显现为具有不同颜色的斑点。此方式也适用于预先涂布专用荧光指示剂F366的色谱板。荧光物质和物体会在明亮的背景中显现为黑色或彩色斑点。F254等其他荧光指示剂,在这种紫外线照射下不发射荧光,因而不会干扰测量。
- · 短波紫外线(UV-C) (254 nm)光照基本上只适用于薄层色谱图。如果色谱板预先涂布F254荧光指示剂,则吸收此种光线的物质在明亮背景下显现为黑斑。

用户可以评估记录下来的板图像,生成报告或导出数据,以便使用其他系统进行进一步处理。

开发TLC Explorer旨在供实验室使用,须由训练有素的用户执行薄层色谱分析,且仅应使用最大尺寸为20 × 20 cm 并且完全干透的TLC色谱板。

请注意,除上述设计用途外的任何其他用途将使保修条款无效。

1.2 安全与处理说明

TLC Explorer的开发,始终以操作安全性为主要目标。仅当严格根据本手册中的说明操作本设备时,才能保证安全。 并且,与本手册说明的安全性简介相冲突的任何操作,都将使保修条款无效。因此,所有负责安装、使用和维护本设 备的人员,都必须在使用TLC Explorer之前仔细阅读本操作手册,并将手册置于手边以便随时查阅。

简介

1

3

9

10

11

版本1.0-2024/02

1

- 5
- 6

7

8

9

10

11

12

1.2.1 在以下情况中,切勿操作TLC Explorer ...

在为TLC Explorer连接电源之前,请确保**杜绝**以下任一情况:

- 设备已经在不良条件下储存了一段时间。
- 设备在运输过程中已经损坏,或经不正确的处理后已经损坏(例如,在运输过程中坠落或接触了液体)。
- 外壳、前面板组件、背板、抽屉或底板的任何部分已经浸湿、损坏或变形。
- · 防紫外线检测窗丢失、损坏或变形。
- · 设备未放置在干燥、平整和稳定的表面上。
- 抽屉无法完全拉出,或无法正确关闭。
- · 未从设备中取出运输过程中防护用的泡沫板。
- 背板的安装不正确,所有四个螺钉都未拧紧。
- · 任何连接组件(USB-C、USB-A、以太网和WiFi天线)或线缆已经浸湿、损坏或变形。
- · 电源不是推荐的两种适配器之一(USB-C电源适配器DELL HA65NM170,或移动电源DELL PW7018LC)。
- · 设备温度明显低于环境温度,此时,空气中的湿气会冷凝并损坏电路(请等候设备达到室温)。
- · 设备所处环境温度不在规定的 15 40°C范围内,或相对湿度未达到20 80%。
- 设备目前或曾经暴露于酸性、碱性、溶剂蒸汽或其他可能导致腐蚀的气体中。
- · 液体或其他物质泼溅到设备表面上或内部。如果发生上述任何一种情况,都请立即移除设备。请确保于再次使用前,设备未损坏且为完全干燥状态。

如果预见到您设备出现上述任何问题,切勿连接、通电或进行操作。有关进一步操作的任何疑问,请联系您的TLC Explorer供应商,或通过图1.1中所示链接访问TLC Explorer服务网站:





1.2.2 安装要求

当您在实验室中安装设备时,请遵守下列指导原则,以保障操作的安全性与可靠性:

- 设备必须在能够支撑10千克重量设备、平整且干燥的实验台上进行安装。
- · 切勿使仪器遭受任何强烈震动或冲击。
- · 保护仪器隔绝过多灰尘。
- ·如果适用,使TLC Explorer、线缆和移动电源与热源(例如,燃气灶、电加热器或烤炉)、溶剂等化学物质或可能导 致本设备受损的其他设备保持安全距离。
 - · 仅使用USB-C电源适配器DELL HA65NM170或移动电源DELL PW7018LC为设备供电,将其连接到本设备背面的USB-C端口。
 - · 根据您实验室的电气设备使用规则,将电源连接至距离本设备2米以内且具有接地保护的插口。
- 用户场所的线电压必须满足电源适配器上指定的规格。
- 切勿将本仪器放置于产生高强度电磁场的设备附近,例如电焊设备、高频熔炉和杆式变压器等。
- · 本仪器连接的电源线须不受突变或电压波动的干扰。
- · 如果电机驱动的设备 (例如, 搅拌机或振动器) 必须与TLC Explorer共用电源线, 请确保安装降噪装置。
- · 确保线缆的布线既不会对线缆操作人员、附近的工作人员或本设备产生风险,也不会导致它们被从插口中意外 拔出。
 - · 设备前方必须留有足够空间,以便能够完全打开抽屉和安全便捷地处理底板。
 - · 设备放置的高度,应使用户能够方便地观看检测窗,并避免外部光源在前面板玻璃上的反射光,以防干扰透过检测窗进行观查。
 - · 设备仅应在满足所有官方实验室安全性要求的环境中使用,且仅应以满足此要求的方式进行操作。

1.2.3 紫外线照射

评估TLC色谱板所使用的用于激发荧光的紫外线照射,会导致眼睛和皮肤受伤。因此,大部分国家/地区的安全性规定,都要求保护眼睛和皮肤免受有害紫外线照射。TLC Explorer内建多种安全措施,能够确保在按照本手册说明使用时不会从设备泄漏长波或短波紫外光。无论何时,不得尝试规避这些安全措施!

- · 仅通过防紫外线检测窗观查紫外线照射下的TLC色谱板。
- ·确保检测窗没有损坏或变形。
- · 在激活紫外线光照之前,请检查抽屉和背板已正确关闭且未损坏。
- · 切勿在TLC Explorer设备外部操作光照装置。
- · 操作紫外光时,请遵守一般实验室安全性规定。

3

4

5

6

9

10

11

12

- 3
- 4
- 5

- 8
- 9
- 10
- 11

12

1.2.4 一般操作注意事项

- 确保在拔掉电源之前,设备至少有30秒的时间彻底关机 这一点与台式机电脑类似。
- · 仅记录无溶剂残留的TLC色谱板 请确保TLC色谱板在上一次色谱展开或衍生化步骤之后已经完全干透。
- · 切勿在底板上放置高于8 mm的物体,这是因为关闭抽屉时可能损坏设备,并且会在图像捕捉时失焦。
- 切勿握着抽屉把手提升或搬运设备。相反,应握住底部的橡胶脚,防止其从您手中滑脱。
- 一旦设备停用,应关闭抽屉以防止损坏和污染。
- ·运输TLC Explorer时,必须关闭TLC Explorer并且移除底板。不应插入任何TLC色谱板。
- · 仅可在USB-A端口上插接FAT32格式的USB棒,进行备份和更新。切勿使用外部硬盘和任何通过线缆连接的其他 介质。
- · 除更换光照装置外(参见第10.4节),仅应由获厂商授权的专业人员打开、调试或维修TLC Explorer。如未遵守这一点,将使所有保修条款失效。
- · 请保留原始包装(包括内衬),以便将来必须运输时保护仪器不受撞击;尤其是在申请保修时,这一点很重要。请 注意,由于不当运输所造成的损坏不在保修之列。

1.2.5 用户管理与信息安全

假定设备放置在实验室中,并且仅限获授权的人员实际接触它。为优先采用平稳高效的实验室工作流程, TLC Explorer不限制特定用户的访问或使用密码或加密技术保护数据。用户账户主要用于将记录的数据结构化。 没有具特殊权限的管理员账户。

实际操作或通过网络连接至设备的任何人皆可

- ・创建新用户
- ·删除现有用户或以现有用户的身份登录
- ·访问和删除所有存储的数据
- ·重启或关闭设备
- · 更改设备设置(包括连接参数)
- ・执行恢复出厂设置。

根据您选择如何在IT环境中集成TLC Explorer (参见第8节),大量用户可以访问您的一台或多台设备。如果 TLC Explorer接入了LAN或WiFi网络,一般来说,能够访问该网络的任何人,都可以打开您设备的用户界面并以现有 用户或新用户身份登录(参见第3.4节)。如果您很重视保密性,请将保密数据复制到您的台式机或平板电脑中,然后 从TLC Explorer中移除相应数据。

- · 计划并仔细测试TLC Explorer的连接设置,确保只有获授权且经过培训的人员具有操作设备的权限。
- · 下载或复制您的数据到存储棒中,然后将其存放在外部储存位置。切勿将TLC Explorer用作长期数据存储介质。
- 仅应用来自厂商经认证的更新。
- · 仅连接您确定不包含恶意软件的USB存储棒(闪存驱动器)。

3

4

5

6

7

1.3 TLC简介

薄层色谱法(TLC)是一种用于定性和定量分析的简单易用和高度灵活的快速分离技术。它非常适合快速识别、筛选和反应监测。其强大的矩阵容错能力和同步分离多个样品的能力,使得TLC具有高时效性和经济性。它广泛用于样品鉴定和定量分析。薄层色谱法几乎可用于分析任何类别的物质,包括农药、类固醇、生物碱、脂质、核苷酸、糖苷、糖类和脂肪酸。

1.3.1 TLC工作原理

常规薄层色谱分析流程简单易用,如图1.2所示。



图1.2:TLC流程:层的制备(色谱板),样品制备,固定相的选择,液体流动相的选择, 样品应用,展开,衍生化,记录

工作流程中的关键步骤是色谱分析过程。薄层色谱法(TLC)遵循经典色谱原理,基于混合物各组分在固定相和液体 流动相之间的亲和力差异进行分离。固定相是惰性载板表面(通常是玻璃、塑料或铝)涂布的薄层吸附材料层(通常 是硅胶或氧化铝)。将样品点样到薄层板(起始线)一端,并垂直插入装有有机溶剂(流动相)的密封室中。流动相通过 毛细作用力在板上向上移动,样品各组分因其对固定相和流动相的亲和力差异而迁移不同的距离。当溶剂到达板的 上三分之一高度时,将薄层板从展开室中取出并干燥。分离组分呈现为板上的斑点,以此计算各组分的比移值(R_f)。

1.3.2 比移值R_f

R_f定义为单个组分移动距离除以溶剂移动总距离;它的值总是在0到1之间。为避免使用小数点,比移值也可以乘以 100,用hR_f表示。

通常化合物与固定相吸附剂的结合越强,它在板上的迁移就越慢。由于TLC吸附剂一般是极性的,因此非极性化合物在薄层板的移动较快,R_f值较高;极性化合物一般移动较慢,R_f值较低。理想的流动相会把所有组分移离基线,最终的R_f值在0.15到0.85之间。

若分离的样品组分是无色的或者说对紫外线照射没有反应,因而不发射荧光,则在进行分析和记录之前,应通过适 合的检测试剂完成分析物的衍生化或可视化。

TLC Explorer记录系统为可靠的TLC色谱板分析、数据处理和数据存储提供便利的解决方案。

1.4 关于本手册

以下各节的编写是为了在了解一般工作流程和软件使用(第4节)之前,首先熟悉设备硬件(第2节)并指导您将 TLC Explorer投入使用(第3节)。请阅读第10.1节,了解如何清洁设备的建议步骤。

您可视需要参考后续章节,它们将提供有关设备设置(第5节)、用户和数据管理(第5节和第6节)以及故障排除 (第9节)的技术细节。

第8节"连接TLC Explorer"旨在帮助IT人士将一台或多台设备集成到您的IT环境中。

如果不确定术语在本手册语境中的具体含义,请参考术语表(第11节)。您还可以在术语表中了解到更多有关用于生成和处理数据的技术与算法的信息。

在涉及按钮或连接器等特定设备元件时,本手册一般会引用在"设备概述"(第2节)中使用的ID。通过查看图2.1、 图2.2和图2.3中的ID,能够轻松在设备上找到对应的元件,如以F开头的ID代表"前面板"上的元件,B开头的ID代表" 背板"上的元件,以P开头的ID代表"板"。

11

2 了解TLC Explorer

2.1 TLC Explorer的正面



图2.1:识别设备正面的元件

2.1.1 F1 - 电源按钮

要启动或关闭设备,请用手指按住电源图标F1(触控按钮)至少两秒钟。在设备启动或关闭过程中,图标的背光将闪烁几秒钟。

暗淡白光指示设备连接到了电源,但尚未关闭。更明亮闪耀的图标则指示设备已开机且为使用准备就绪。 请牢记,要关闭设备,须按下电源按钮(或在软件中选择Shutdown(关机))并等待图标停止闪烁,然后拔下电源 插头。

2

3

1

5

6

8

9

请注意,通过上下滑动滑块F8,即可通过检测窗F7观查被照射的色谱板。光照按钮仅用于为目视检查而打开/关闭光源,不用于捕捉图像。

将手指放在按钮标签"Vis"、"366"和"254"开关右侧的圆圈上,可打开或关闭对应的光照。但一次只能触发一种光

如果紫外光可能从设备中泄漏,则安全功能会自动关闭光源并禁用这些按钮,例如,未正确关闭抽屉F5或未正确安 装背面罩板B7。

Vis: 按钮F2,标为"Vis",开启白光LED光照。

2.1.2

F2, F3, F4 - 光照按钮和指示灯

366: 按下F3, 切换为长波紫外线LED光照, 可查看以366 nm波长照射时显示荧光的物质。

源。圆圈后面的背光,指示光照已经打开-用户应按下按钮或通过软件控制来手动捕捉图像。

254: 按下F4,切换为254 nm波长的短波紫外线LED照射。使用此光源可查看吸收这一波长的物质,它们在预先涂 布F254荧光指示剂的TLC色谱板上显现为黑斑。

2.1.3 F5, F6 - 抽屉和抽屉打开指示灯

轻按抽屉把手的中间部分以打开抽屉F5。如果设备已开机,您会看到电源按钮F1下方的图标F6亮起红光。在任何时候,当传感器指示抽屉未正确关闭,即会亮起红光。抽屉打开后,无法打开光源和拍摄图像。 要在底板上放置TLC色谱板或从底板上取下色谱板,需要完全打开抽屉。

请将多个色谱板整齐排列在底板上,然后轻按抽屉把手的中间部分直至您感觉到轻微的阻力。抽屉最后几厘米的滑 动将由抽屉自主完成关闭。专门的阻尼系统能够让您正确关闭抽屉而不会移动色谱板。

在系统记录图像时打开抽屉,会导致软件显示错误信息并退出捕捉过程。

2.1.4 F7, F8 - 检测窗和滑块

打开所需光源(F2、F3或F4),然后向下轻按滑块F8,透过检测窗F7观查每个TLC色谱板,是在紫外光或可见光照射下查看TLC色谱板的快速安全方式。

滑块仅能打开大约半个检测窗大小的区域。这一设计减少了用户通过窗口进行观察时入射杂散光的干扰。通过上下 滑动滑块,便可查看底板的不同部分。

版本1.0-2024/02

1

2

3

4

5

6

7

2 了解TLC Explorer - 2.2 TLC Explorer的背面

2

3

4



图2.2:识别设备背面的元件

2.2.1 B1 – 用于插接电源的USB-C端口

背面右上角的USB-C端口B1仅用于连接电源,不用于数据传输。在连接任一推荐的电源后(USB-C电源适配器DELL HA65NM170 A2或移动电源DELL PW7018LC A4),前面板上的电源图标F1应呈暗淡白光亮起,指示设备已为开机做好准备。

2.2.2 B2, B3 - USB-A端口

您可使用左侧(B2)或右侧(B3)的USB-A端口连接存储棒(闪存驱动器)来备份或更新软件。切勿使用外部硬盘和任何 通过线缆连接的其他介质。请确保仅使用无恶意软件的存储棒。

2.2.3 B4 - 以太网端口

根据您的喜好,您的TLC Explorer可通过WiFi(主要是指平板电脑和移动设备)或以太网(例如,局域网/LAN中的台 式机电脑)连接至外部设备。要使用以太网选项,须将常规5类线或6类线连接至RJ45端口B4并视需要按照第8.2节 中的建议配置TLC Explorer和您的局域网。

9

11

3

4

5

7

12

2.2.4 B5 - WiFi天线连接器

每台TLC Explorer均出厂配备一个单独的WiFi天线A1。为将天线连接至设备,请将天线基座上的接头对准背板左上 角的RP-SMA连接器B5,然后顺时针旋入拧紧。您可调整天线来优化信号接收和占用的空间。 有关如何配置WiFi连接的详细信息,请参见第8.1节或第8.3节。



图2.3:安装WiFi天线

2.2.5 B6-序号

背板上的贴纸B6(在USB-C端口的旁边),用于通过其标识的序列号识别特定设备。它应与软件里Network(网络)下Settings(设置)页面中显示的序号一致。

序号下方的二维码包括连接至完整设备标签的链接,完整设备标签通常位于设备的底部。

2.2.6 B7 - 背面罩板

在TLC Explorer的背部,是由4颗花纹螺钉紧固背面罩板B7的。该罩板仅可在清洁设备内部(参见第10.1节)或更换 光照装置(参见第10.4节)时打开。紧固罩板的全部4颗螺钉,以便拆装。 请注意,如果传感器检测到背板未正确安装,则安全机制会禁用所有光源,因而将无法获取图像。

2.2.7 B8 - 光照装置

光照装置B8仅在拆除背板B7后可见(参见图10.3)。其内装不同光照类型的LED光源,是设备的唯一可更换部件。 请检查您的光照装置是否需要更换,有关如何安全地更换,请参见本手册第10.4节。

3

4

5

2.3 TLC Explorer的底板



图2.4:识别底板上的和与底板有关的元件

2.3.1 P1 - 底板区域

将您计划直接记录的TLC色谱板放置在底板区域P1上,色谱板下面不需要纸张或箔纸等任何其他材料 – 自动检测 色谱板所需的黑色涂层是光学仪器的一部分。请确保色谱板平稳放置在底板区域上,切勿伸出底板边框。切勿在底 板区域上放置高于底板边框P2(通常为8 mm)的物体,这样做在关闭抽屉时可能损坏设备,并且会导致图像捕捉时 失焦。

底板区域足以容纳一个20 × 20厘米大小的TLC色谱板或是多个较小的色谱板。要使设备能够自动检测多个色谱板,放置它们时须相互间隔几毫米,以使设备能够检测到在每个色谱板周围都有至少2毫米宽的黑色涂层。 根据喜好,您可以在底板位于抽屉中时,将TLC色谱板放置在底板区域上,或者取下底板放到别的地方,再将色谱板放置在底板区域上,然后将底板连同布置好的TLC色谱板插回设备中。

底板区域四周的隐藏式抓手,使用户能够在不碰触TLC板涂层表面的情况下,从底板上提起多个TLC板。

2.3.2 P2 - 边框

底板区域P1的四周环绕着边框P2,大为方便了底板的处理和运输。在将底板插入抽屉时,不应在底板区域上放置高于边框高度的物体。

2.3.3 P3, P4 - 对准销和底座

底板区域下方的对准销P3,有利于用户在抽屉上正确定向和对准底板。关闭抽屉之前,请确保对准销P3已经插入抽屉底座P4中,以便底板平稳就位在抽屉中。

9

10

11

3

4

5

6

7

8

2.4 附件

2.4.1 A1 - WiFi天线

为提供可靠的WiFi连接,TLC Explorer出厂配备了一个外部天线A1,用户需要将其安装在连接器B5中(有关详细信息,请参见第2.2.4节)。

2.4.2 A2, A3 - 电源适配器和电源线



图2.5:电源适配器和电源线

设备包装中包括一个USB-C电源适配器A2和一套用于将适配器接入电网的电源线A3。请选择适合您本地电源插口的电源线。该电源适配器的输入交流电压为100到240 V,频率为50或60赫兹。

在将电源适配器接入电网之前,请检查所选电源线是否不会遭受突变或电压波动的影响(另请参见第1.2.2节)。一旦 电源适配器接入电网,请将USB-C连接器插入设备背面的插座B1中。如果前面板上的电源指示灯F1亮起,则说明设 备已准备好启动。



图2.6:移动电源 - 不包括在设备交付清单中

移动电源A4不在出厂交付设备清单中,但是可以单独购买(参见第10.3节),它能在没有可用市电插口时为 TLC Explorer供电。充满电后的移动电源可为TLC Explorer提供充足电力,保持续航约30个小时。 使用移动电源而独立于电网时,用户将获得更高的便携性,应始终以对待光学测量仪器的态度给予设备适当照护。 有关详细信息,请参见第1.2节。

版本1.0 - 2024/02

3.1 拆除包装与交付清单

TLC Explorer的制造和检测符合电子仪器的相关指导原则及规范(有关详细信息,请参见第12节),出厂时经过安全测试,处于技术完好状态。如果发现任何可能危及设备的包装损坏,请记录相关情况并联系您的供应商。如果您怀疑符合任何第1.2.1节所列出的标准,切勿操作TLC Explorer。



图3.1:包装概览

强烈建议保留包装材料以便日后运输和/或储存设备,尤其在申请保修时这一点很重要。 打开纸板箱后,首先请取出包括底板在内的所有附件和顶部纸质防护物。确保在取出设备前,取出所有零散的部件。 转动纸板箱,让玻璃前面板正对您,准备好从箱中拿出设备后放置它的适合空间。从相对的两个方向握住设备底部, 然后提起重量约为10千克的TLC Explorer。移除所有运输固定物(如固定抽屉的胶带)和防护箔纸。

检查交付的货物包括下列项目:

- · 一台TLC Explorer,其内装光照装置(检查光照装置,参见下文第3.3节)
- · 一个底板
- · 一个WiFi天线
- · 一个USB-C电源适配器,及适合不同插口类型的八条电源线
- · 一个防尘罩
- 两张贴纸,包含两个二维码 用于更方便地连接设备(参见图3.2)
- 一本名为《TLC Explorer Documentation System Getting started 》(TLC Explorer记录系统 入门指南)的手 册和一份安全附录

11

3.2 适合安全操作的设置

在为您的设备选择适合的安装地点之前,请仔细阅读第1.2.2节中的安装要求。检查插入了底板的抽屉F5是否能够 完全打开,所有计划内用户是否能够以安全、便捷和符合人体工学的方式透过检测窗F7观查TLC色谱板。尤其需要 考虑不同的人体高度和设备玻璃前面板的干扰性反射光。

请将一份用户操作指南放在TLC Explorer的附近,以便随时查看,例如您计划用于控制设备的平板电脑或台式机电脑中的指南电子版本。

3.3 插电和'首次查看'底板和色谱板

挑选适合电源插口的电源线A3,将其连接到电源适配器A2和电网,然后将USB-C连接器插入端口B1中。您应在设备 前面板上看到电源图标F1亮起。用手指轻按电源图标约两秒钟,直至图标开始闪烁,这指示设备正在启动。约30秒后 图标停止闪烁并变得更明亮,此时,设备已为操作准备就绪。

打开抽屉F5,插入底板。具有宽外缘和压花标志的一侧应朝外,即,该侧应与抽屉把手毗邻。检查对准销P3是否已经 插入抽屉底座P4中,以及底板是否平稳就位在抽屉中。要关闭抽屉,应轻按把手的中间部分直至您感觉到轻微的阻 力。抽屉最后几厘米的滑动将由抽屉自主关闭。专门的阻尼系统能够让您正确关闭抽屉而不会移动色谱板。

一旦红色的抽屉打开指示灯F6熄灭,用手指短暂按住标签 'Vis' F2旁边的圆圈,打开内部的白光光源。向下滑动滑块F8,透过检测窗F7看向里面,您应看到被照射的底板。

关闭光源(触按F2),打开抽屉并将一个TLC色谱板放置在底板区域P1上。再次关闭抽屉,透过检测窗观查不同类型 光源照射下的色谱板(触按F2、F3或F4)。

如果您的TLC Explorer运行表现异于上文描述,请在操作设备之前参考第9节来识别和解决问题。

1

4

9

10

11

2

3

4

5

6

7

8

3.4 连接笔记本电脑

将多种用户设备(笔记本电脑、台式机电脑、平板电脑或智能手机)连接到TLC Explorer的方法各异,请根据第8节进行配置。推荐的基本方法是将笔记本电脑连接到TLC Explorer自己的WiFi网络。此方法的设置很快捷,只需点击几下鼠标,即可将生成的图像和数据直接下载到您正在使用的设备上。



图3.2:通过扫描设备标签上的这两个二维码就能轻松连接

在笔记本电脑(或具备WiFi的台式机电脑)上打开网络设置,搜索以'tlc-explorer'开头且后跟序号B6的网络名称 (SSID)。选择'WPA/WPA2'为安全性设置,输入步骤1标签(参见图3.2)上显示的密码,建立连接。连接平板电脑或智 能手机的方法则更加简单,只需扫描设备提供的贴纸上的二维码1。您可能会看到提示所选WiFi没有互联网连接的 警告,根据设计这属于正常情况。

一旦建立了WiFi连接,就可以在您的网页浏览器中打开用户界面。尽管可以使用所有通用、最新款的浏览器,但推荐您使用Chrome或Firefox浏览器,以获得最佳的用户体验。

要打开TLC Explorer用户界面,您可以扫描贴纸上的二维码2,或在浏览器地址栏中输入"http://tlc",或用设备名代 替 'tlc' (有关详细信息,请参见第8.2节)。

如果浏览器难以搜索到"http://tlc"(例如,由于设备通过移动数据网络连接到互联网),您还可以输

入"http://192.168.200.1"。有些浏览器会强行建立加密连接,可能会自动为输入的地址加上"https://"前缀,然后 会生成提示网站可能不安全的警告信息。浏览器无法知道您正在访问受保护局域网中的TLC Explorer,其风险远比 从互联网打开网站小得多。可以忽略警告继续操作,也可以在浏览器设置中禁用https-only(仅https),然后在浏览 器地址栏中再次手动输入"http://tlc"。



图3.3:初次打开的TLC Explorer用户界面

建立书签和/或从浏览器菜单(通常位于浏览器窗口的右上角,标记为三个竖向的圆点)中选择菜单项Add to Home Screen (添加到主页屏幕),这可以显著简化未来打开某网站的操作。

3.5 创建首位用户

通过选择语言和输入首位用户名,开始个性化设置您的TLC Explorer。由于用户名用于内部命名文件和目录,所以不得使用任何特殊字符或符号。如果输入了无效字符,则栏位下方的线将变为红色,并将在下方显示警告信息。您可以在下一个栏位输入全名,此项非必填项。例如,它可用于识别报告作者用户。

Add a nov			
Add a flev	v usei		
Language			
English		~	
Username			
EgonStahl			
Full name (optional)			
Egon Peter Gus	tav Stahl		
 Device 	and network settings		
 ✓ Legal te 	erms		
Accept legal	terms		
The device is	used in the USA or Canada		and the second se
	Add user		

图3.4:添加首位用户

目前可保持Device and Network Settings (设备和网络设置)为默认值。日后可随时在Settings (设置)页面进行修改。

在勾选Accept Legal Terms (接受法律条款) 左侧的复选框之前,请展开和仔细阅读法律条款。操作设备和软件需要您同意这些条款。

如果您的设备位于美国或加拿大,需要您勾选相应的方框。

点击Add User(添加用户)后,即完成了初始设置,如此设备就已为使用准备就绪。

2

版本1.0 - 2024/02

您将被自动转至如图3.5所示的标准欢迎屏幕。选择您创建的用户,并点击Login(登录),便可开始捕捉TLC色谱板的图像。

TLC EXPLORER DOCUMENTATION SYSTEM Supelco [®] Analytical Products	Welcome!	≡
	Username	
	EgonStahl	
		Login
© 2023 Merck KGaA, Darmstadt, Germany and/or its affiliates. All rights reserved.		

图3.5:欢迎屏幕,其提供用户的选择

3

4

5

6

7

8

9

10

关闭设备

3.6

和所有台式机电脑一样,在拔下电源插头之前,需要正确关闭TLC Explorer。否则,可能损坏设备。

为正确关闭设备,请点击主菜单图标(右上角的三条横线),从显示的菜单中选择Shutdown(关机)。在如图3.6所示的对话框中,可以选择关闭设备、重启(即,重新引导)设备或仅重启软件应用程序。您也可以触按电源图标F1约2秒钟,但需要在用户界面中确认关机(在所有活动的用户会话中都将自动显示前述的对话框)。

如果选择重启设备或应用程序,只需等待约1分钟,直至图3.5所示的欢迎屏幕重新显示,这时设备已经再次准备就 绪。

 Warning X	_
Are you sure you want to shutdown the system? Unsaved data will be lost. Please make sure all data is saved before proceeding.	
Cancel Software restart Device restart	
Shutdown	

图3.6:关闭或重启设备

目前可保持Device and Network Settings (设备和网络设置)为默认值。日后可随时在Settings (设置)页面进行 修改。

在勾选Accept Legal Terms (接受法律条款) 左侧的复选框之前,请展开和仔细阅读法律条款。操作设备和软件需要您同意这些条款。

如果您的设备位于美国或加拿大,需要您勾选相应的方框。

点击Add User(添加用户)后,即完成了初始设置,如此设备就已为使用准备就绪。

12

4 操作指南

4.1 典型工作流程

无论是仅快速一览TLC色谱板、捕捉和导出色谱板图像(可能带有注释)进行记录,或需要使用完整报告进行定量分析,TLC Explorer允许您按需操作。

因此,不需要遵守一成不变的流程。您可选用多种不同的工具并组合使用这些工具来视需要处理数据,例如,为了调查不同的研究问题,您甚至可以用多种方式评估同一个TLC色谱板。

下图4.2概括了不同的记录和评估步骤。有关这些步骤的详细信息,请参见本节中的相应部分(按文字左侧的索引编 号查阅对应章节即可)。

在图4.2的右侧,您将看到带有Image(图像)、Plate(板)和Evaluation(评估)字样的竖条。这与软件用户界面(参见 图4.1)顶部的进度条相对应,本手册用了3页的篇幅介绍处理步骤:

- · Picture (图像): 捕捉新图像或加载现有图像集。
- · Plate(板): 命名一个图像集中的每个色谱板、添加信息和定义轨迹。
- · Evaluation(评估): 注解每个色谱板图像、绘制和比较轨迹密度图或执行定量分析。



图4.1:主要步骤的进度条和右上角的主菜单(在竖向3个横条上点击)

3

4

5

6

8

9

10

11

4 操作指南 - 4.1 典型工作流程



4.1.1 主菜单

点击右上角的3个横条,打开主菜单。通过主菜单,可访问在工作流程图中未占据一席之地的功能。除只提供 Settings(设置)和Shutdown(关机)功能的欢迎屏幕外,主菜单浓缩了图4.3显示的所有页面和选项。有关详细信息,请查阅这些选项右侧列出的章节。



图4.3:包含本手册相关章节的主菜单

4.2 准备设备

在将TLC Explorer接入电网或移动电源后,电源图标F1后面透出的暗淡白光表示设备已准备好启动。用手指轻按图标约2秒钟,等待设备准备30秒左右,当图标开始闪烁,就说明设备已为操作准备就绪。 把一个TLC色谱板放置在底板上,关闭抽屉F5。当红色的抽屉打开指示灯F6熄灭,您将看到'Vis'F2旁边的圆圈呈现短暂的白光闪烁 – 在此期间,系统完成快照并在软件中显示色谱板的预览。

2

9

10

4.3 目视检查

前提条件:

- · TLC阅读器打开;电源图标F1以亮光亮起(不是闪烁或暗淡白光)。
- · 抽屉正确关闭,红色的抽屉打开指示灯F6熄灭。
- · 没有用户正在同时捕捉图像(另请参见第4.4节)。

选择您要使用的光照类型,请用手指短暂按住相应的按钮(F2、F3或F4),直至圆圈亮起。然后轻按滑块F8,打开检测 窗F7。根据观查角度,您可上下滑动滑块,观查TLC色谱板的相应部分。 通过触按相应按钮,还可切换至其他光照类型,也可通过再次按下同一个按钮关闭光照。

4.4 捕捉图像

前提条件:

- · TLC阅读器打开;电源图标F1以亮光亮起(非闪烁或暗淡白光)。
- · 抽屉和检测窗正确关闭,红色的抽屉打开指示灯F6熄灭。
- ·最小尺寸为2.5 × 7.5厘米的多个或最大尺寸为20 × 20厘米的单个TLC色谱板平稳就位于P1底板区域上。
- · 同一时间仅一位用户可以捕捉图像。尽管其他用户也可加载和处理图像集,但若同时捕捉图像将会导致提示错 误信息。

假定您已经按照第3.4节中的建议,连接了平板电脑或智能手机,则只需点击桌面图标或在浏览器中保存的书签, 即可打开如图3.5所示的欢迎屏幕。选择您的用户名,并点击Login(登录)。

4

5

6

7

8

9

10

11

在每次登录(如图4.4所示)后打开的页面中,可以捕捉新图像,或者加载现有图像集(参见第4.5节)。在屏幕左侧 (如屏幕尺寸较小,则在顶部),您将看到在关闭抽屉时捕捉的低分辨率底板图像。只需点击该图像,即可刷新图像。



图4.4:捕捉图像或加载现有图像集

从下拉菜单中选择Capture new Image Set (捕捉新图像集)选项,并在文本栏位中输入新图像集名称。此名称将用 作图像文件名的一部分,因此不是任何字符都被接受的。而且,此名称需要与该用户保存的其他图像集名称有所不同。如果图像集名称无效,软件会提示您原因。

复选框让您可以选择图像捕捉所用的光照类型。请注意,根据默认设置会激活 'VIS' (白光);待选中其他光照类型 后,将取消选择VIS。

展开Use manual Exposure(使用手动曝光),您可以为每种选定的光照类型设置曝光时间,单位为毫秒(参见 图4.5)。点击相机图标,系统将按照所选曝光时间拍摄低分辨率图像,并显示此图像作为预览。点击Auto(自动)按 钮,可要求系统为色谱板拍摄一系列图像,并自动决定最佳曝光时间。请知悉,拍摄过程可能耗时几秒钟,当您看到 预览图像更新后,才可继续操作。

点击Capture Now (立即捕捉), 仪器将开始捕捉图像。通过进度条上方的文字和设备上亮起的相应光照按钮 (F2 - F4), 您可以知道当前正在拍摄哪些图像。 所有光照类型生成的图像,将一起保存为一个Image Set (图像集)。然后,系统会在多个图像中搜索色谱板。一旦检测到所有色谱板、旋转了所有色谱板 (将其底部边缘水平对齐)并提取为多幅单独的图像进行进一步处理后,用户将 被自动转至下一步进行板编辑 (参见图4.6)。

如果仅希望使用捕捉的图像(或提取的色谱板图像),您可以通过文件管理器下载刚刚拍摄的图像,从而忽略所有进一步的处理(有关详细信息,请参见第4.14.1节)。还有一种可选方法,点击Plate(板)页面底部的Download image (下载图像)按钮,保存当前在预览区显示的图像。未来任何时间均可加载和处理图像集。



图4.5:为选定的光照类型手动设置曝光时间

4.5 加载现有图像集

前提条件:

・您已捕捉、导入或复制了图像集(参见第4.14.2节)。

要打开和处理此前保存的图像集,请从Image source (图像来源)下拉菜单中选择Load Image Set (加载图像集), 然后点击 – 最初是空白的 – Select an Image Set (选择图像集)下拉菜单。 显示下拉菜单,它允许您逐一预览图像集中的图像,或是预览整个图像集(如图4.6所示)。

点击Use Image Set (使用图像集)来确认您的选择。系统将自动转至下一步进行板编辑(参见第4.6节)。 用户只能打开自己的图像集。要使用属于其他用户的图像集,您需要在文件管理器中创建自己的拷贝(参见 第4.14.2节)。



图4.6:预览和打开现有图像集

4.6 选择和编辑色谱板

前提条件:

- · 您已捕捉了图像或加载了一个图像集。
- · 至少找到了一个板。

在评估轨迹之前,您可以分别命名和编辑在Image Set (图像集)中找到的每个板。也可选择添加有关板的信息和修改轨迹属性,构建供进一步分析的基础。

在左侧的预览区(参见图4.7)(或者屏幕尺寸较小时:预览区位于顶部),板已经对齐和被裁切(具有较小的图像页边 空白),并且系统将自动在检测到的轨迹四周绘制边框。

如果您将多个板放置在底板上,点击Show Overview (显示一览),可查看所有板 (带索引编号),然后从顶部下拉菜 单中选择要编辑的那个板。特别是当一个图像集中有多个板时,建议您使用Rename (更名) 输入栏位为每个板命名 一个独特的名称。

2

3

9

11

ОК

ОК

2

3

4

5

6

7

8

9

10

11

12



图4.8:选择光照和频道

· Original (原图),是指RGB彩色图像,

('UV366')。使用此图像,您可以选择特定的频道,即

· Intensity(强度),图像灰度化,

- · Red (红光),仅红光频道的强度,
- · Green(绿光),仅绿光频道的强度,
- · Blue(蓝光),仅蓝光频道的强度。

为改进带伪彩色的低对比度单频道视图的可视性,可激活Channel(频道)下拉菜单右侧的Heatmap(热图)复选框。

您可采用+90°和-90°按钮来旋转板的图像。旋转板图像还将删除所有轨迹,并由在新方向上自动检测到的轨迹取代

原来的轨迹。使用此功能,可重置轨迹搜索,例如,板的轨迹宽度与默认值不同。如果需要,展开Tracks(轨迹)区,设

ок

图4.7:选择要编辑的板

在板旋转按钮的下方,是两个分别名为Illumination(光照)和Channel(频道)的下拉菜单,在其他处理步骤中也会 遇到它们。设备会分别针对捕捉图像而选定的每个光照类型(参见第4.4节)创建一幅8位RGB图像(即,它保存了红、 绿、蓝光的强度,是彩色图像)。Illumination(光照)使您能够选择特定的板图像,例如,白光('Vis')或长波紫外线

置适合的轨迹宽度,然后来回旋转板,这会发起新的自动轨迹检测。

 \odot

为方便日后参考,在捕捉图像之前,可视需要在Info(信息)区输入有关板及其制备的可选详细信息,以利生成报告。 这包括分析所使用的板类型、溶剂和染色相关信息,以及您希望与板数据一同存储的说明和评论。

也可编辑板图像以去噪或增强对比度(参见图4.9)。为防止原始数据被覆写,需将修改后的图像单独保存,方法是在 New processing name(新处理名称)下方输入名称并点击其右侧的按钮。这样一来,在所有未来处理步骤中,均可 从Illumination(光照)下拉菜单中选择修改后的图像。请注意,Quick clean(快速清洁)和Sensitive clean(敏感清 洁)算法,能够去除长波紫外线光照(波长366 nm)所捕捉图像中的杂质和污染,也可能会生成伪影。

M		TLC EXPLORER		08:40 AM 🗮
	Picture	DyeMix lot 42	Evaluation	
and the second		A start for the start	1: DyeMix 20x10 cm	~
			Rename	
			DyeMix 20x10 cm	
-			Show overview Correct plate orientation +90° Illumination Channel UV366 VOrigina	-90°
			∽ Info	
			 Post processing 	
			Denoise Quick clean	Sensitive clean
			Enhance contrast	
			20	255
			New processing name	200
			A suffix for the new image file	
	_			
	图4	4.9:图像后期处埋,溶剂前沿可见的示(例	

3

4

8

- 3
- 4
- _
- 5

12

4.7 定义轨迹

前提条件:

- · 您位于Plate(板)页面(参见顶部的进度条)
- 板的方向正确(旋转板将重置您对轨迹所做的更改)。

此步骤允许您添加、删除、移动和命名轨迹。轨迹的定义将用于在下一步开始的新的评估。如有必要,您日后可以回 来编辑轨迹,创建带有修改后轨迹属性的新的评估。但为确保数据完好性,任何对轨迹的更改都不得影响现有的评 估。

自动轨迹检测会考虑到所有捕捉的图像 – 也是为了检测仅在紫外线光照下可见的物质 – 然而,可能不会总是发现 所有轨迹都是正确的。如果发现默认轨迹宽度与板的轨迹宽度不匹配,请调节Track Width (轨迹宽度) 滑块,然后来 回旋转板,从而发起新的自动轨迹检测。您可轻松删除错误检测的轨迹,只需点击特定轨迹行上的回收站图标即可。 添加轨迹的方式有两种:在点击Add Track (添加轨迹) 后,您可以

- · 输入距离左板边缘的轨迹位置,单位为毫米,并点击Done(完成)进行确认,或
- · 直接点击预览图像中的轨迹。

在上述两种操作中,添加完所有需要的轨迹后,都需要再次点击Done(完成)进行确认。



图4.10:添加和删除轨迹

您也可以通过更改Center(中心)值(同样,是距离左板边缘的位置,单位为毫米),来移动每个轨迹并为轨迹命名一个独特的名称。

为获得最大的灵活性,可为所有轨迹共同设置基线(起始线)和溶剂前沿,然后为每个轨迹单独微调。一般来说,可以 使用几个滑块来优化调节居中的轨迹基线和溶剂前沿位置(参见图4.11)。假定溶剂前沿在板边缘上的位置较低,可 通过点击相应行左侧的箭头微调第一个轨迹。当取消选择Solvent front(溶剂前沿)前面的复选框后,可以设置该轨 迹的相应值。要再次应用该轨迹的溶剂前沿,只需再次勾选复选框。

点击箭头可打开一览表从而展开轨迹视图(也可以关闭一览表),还允许您为每个轨迹添加说明或评论,记录应用的 样品数量,把轨迹标记为参考(带给定浓度),以便日后进行定量分析。

根据首选的工作流程,日后您还可以为选定的评估来更改基线、溶剂前沿、样品数量和参考属性。

点击Back(返回)按钮,您将退回至捕捉或加载图像集的Picture(图像)页面。请注意,这不会撤销或反转您在Plate (板)页面中所做的设置。点击OK(确定)按钮将带您前往Evaluation(评估)页面(所有与评估相关的轨迹都应预先 定义好)。



图4.11:调节所有轨迹的基线和溶剂前沿,并微调单个轨迹

3

9

11
12

4.8 创建评估

前提条件:

- · 您位于Evaluation(评估)页面。
- 在创建新的评估之前,请检查该评估所使用的所有板的全部相关轨迹是否都已定义完毕。

无论何时创建新的评估后,系统将把板编辑区(参见第4.6和4.7节)中的所有当前板图像和设置复制到另外一个目录中。如果点击了Create new Evaluation (创建新评估)按钮,则日后的修改将仅影响那些在修改后创建的评估(参见 图4.2)。



图4.12:创建新评估

点击板区域底部的OK(确定)按钮,将自动创建名为 'Evaluation_1'(评估_1)的评估。与板相同,您可使用 Rename(更名)文本栏位为其命名一个独特的名称。

另外,也可基于现有评估来创建新的评估。在选择要从此区域顶部下拉菜单中复制的评估后,只需点击 Duplicate current Evaluation (复制当前评估)。对原评估和复制的评估之一进行更改,不会相互影响。 在评估过程中的任何时间,您都可以视需要配置预览区(参见图4.13)。您可以选择希望查看的某个板、该板图像 ('光照')及该图像的特定频道(有关详细信息,参见第4.6节)。而且您可以激活或停用以下项目的显示

・溶剂前沿,

・基线,

- ・ 注释(参见第4.9节),
- · Rf量尺,从基线的0值,到溶剂前沿的1或100(100是因为选择了'使用hRf值'),
- · 一个环绕板的厘米量尺(参见图4.13中的示例)及
- · 环绕所有轨迹区域的定界框。

利用两个滑块可以调节激活的评估的基线和溶剂前沿位置。此区域底部的Capture/Load Image(捕捉/加载图像)和Plate(板)按钮,让您能够导航回至相应页面(可选方法是,点击页面顶部的进度条)。



图4.13:配置用于评估工作的预览区

4

2

1

11

4.9 注解板图像

前提条件:

- · 您位于Evaluation(评估)页面。
- 注释是评估的一部分。如果希望日后编辑注释,请确保已经从此区域顶部的下拉菜单中选择了相应的评估。

您可以在板上标记多个点并添加评论。点击Evaluation(评估)区中的Annotations(注释), 再点击+Add Annotation(+添加注释)以激活此功能。然后在板预览区中点击要添加评论的位置。



在上方表格Annotations(注释)标签下,您将看到与您添加的每个点相对应的各个注释行(参见图4.14)。行的索引编号与您添加它们的顺序相对应。您可以轻松更改这个顺序,通过触按(或鼠标左键点击)并按住某一行,即可在表格中上移或下移该行。在同一表格中,还可以向注释添加文本或从注释删除文本。

虽然在预览中不显示注释文本,但通过按下Export(导出)按钮可以将这些文本写入板图像中它们对应的点旁边,然 后可在点击Download Export(下载导出)之前,选择对话框中的Export images as displayed(导出显示的图像)。 在下载的zip文件中,所有带注释的图像都标记有'_overlay.jpg'后缀。

11

也许您希望日后使用其他仪器来分析相关的板位置,例如进行质谱分析。对于此类应用,注释位置还会列在距离板左下角几毫米的地方(参见Position(位置)列标题下方的下拉菜单)。还可将注释导出为CSV文件(标记有'_Annotations.csv'后缀),这使得在其他设备上输入选定的板位置变得简单。如需导出图像和数据的更多信息,请参见第4.12节。

您将在刚才讨论的上方表格下方看到第二个表格(下方表格)。第二个表格列出了作为Comparison(比较)评估步骤 一部分的斑点(参见第4.10节),此处列出仅作为信息显示。为在预览中区分斑点标记和用户注释,请查找":"-斑点 标记的格式为<track index>:<spot index>,而可自由布放的注释仅显示为一个数字(它们的索引编号)。

4.10 绘制和比较轨迹密度图

前提条件:

- ・ 您位于Evaluation (评估) 页面。
- 在创建评估之前,已在板页面中定义了所有相关的轨迹(参见第4.7节)。

Comparison(比较)标签页下方的工具,允许您绘制和比较轨迹密度图。所选板图像的某一特定频道的密度图(有关详细信息,请参见第4.6节),被视为轨迹宽度上多个密度像素值的平均值或中位数。您可以视需要微调参数,如平滑或线条颜色,还可为斑点添加注释。然后将密度图保存为图像并将此图像纳入文档记录中,或者导出密度图数据以便使用其他工具进行进一步处理。

4.10.1 绘制密度图

为图解Comparison(比较)下方可用的关键功能,请阅读以下逐步操作的示例:

- 1) 从Create new Evaluation (创建新评估) 按钮下方的下拉菜单中选择一个板。
- 2) 点击Comparison(比较)标签页。
- 3) 使用滑块调节轨迹宽度。
- 4) 点击上方表格末尾的+ Add Track (+ 添加轨迹) 按钮。
- 5) 通过点击板图像中的轨迹可以选择一个轨迹,或者从上方表格中的Select a track (选择轨迹)下拉菜单中选择 一个轨迹。
- 6) 在上方表格中的Select an Illumination (选择光照) 下拉菜单中,选择一个用于计算的板图像。请注意,此选择 独立于页面上方的Illumination (光照) 下拉菜单。光照下拉菜单用于设置预览图像。
- 7) 选择图像频道以便计算密度图。选择Original (原图) 会采用与Intensity (强度) 相同的计算基础,比方说,将 RGB值转换为灰度值。然而,密度图下方的轨迹是彩色的,而非灰度图。
- 8) 在板图像上方点击Densitogram (密度图)标签页以查看绘制的轨迹。在板图像下方,您可以看到垂直旋转的轨迹图像,此图像已经过缩放调节,能够显示在密度图的X轴上方(参见图4.16)。
- 9) 为缩放到密度图中,请按下鼠标左键选择欲放大的区域,或是用手指在轨迹图像上方滑动选取欲放大的区域。 下方的轨迹图像将自动适应为显示所选的区域。在密度图上点击一下,可添加一个标尺并显示此R_f值下所有轨 迹的精确值,双击会重置密度图,您又会看到完整的轨迹。

5

8

9

10

11

7

10) 展开Options(选项)区,您能够使用其他控件(参见图4.16),这些控件将应用于此Comparison(比较)的所有 轨迹:

- 将像素行的平均值或中位数值用作密度图值
- 平滑绘制使用的数据,设置过滤器筛选的范围
- · 自动检测和提取基线,即,没有斑点的板的背景级别(通过每个轨迹左侧和右侧的空白区域在本地进行估算)。
- 11) 点击密度图右上角的相机图标,将显示的密度图保存为图像文件。

按照上述方法,您可以添加用于比较的其他轨迹。通过在上述示例步骤1中选择不同的板,甚至可以比较来自不同板 的轨迹(请知悉此种比较的局限性)。为了更好地区分多个密度图,可点击上方表格第一列的线条编辑器,从而选择 线条的颜色和格式。

还可通过线条编辑器右侧的文本栏位来为轨迹更名。请注意,这改变了在所有激活的评估中该轨迹的名称。

4.10.2 添加斑点的名称或注释

利用Comparison(比较)标签页中下方的第二个表格,可以命名(或注解)沿着轨迹和相应密度图的任何位置。激活+ Add Spot(+添加斑点)按钮后,只需点击预览图像中的斑点,或从下拉菜单手动选择一个轨迹,再编辑表格中的hR_f 值(参见图4.17)即可。在相应行的文本栏位中输入斑点名称或评论即可。

然后,此表格中列出的斑点,被视作预览图像和密度图(如果存在特定轨迹的密度图)中的注释。如同常规注释,导出 它们有利于后续处理步骤,像是对单个斑点进行质谱分析。

Parte Parte <th< th=""><th></th><th>09:42 AM 🗮</th><th></th><th>XPLORER</th><th>09:43 A</th><th>м≡</th></th<>		09:42 AM 🗮		XPLORER	09:43 A	м≡
Plate image Densitogram Annotations Comparison Annotations Comparison Quantification Options V Options V Options Task value(well) 0 0 0 0 + Add track 0	Picture DyeMixes	Evaluation	Picture	DyeMtxes	Evaluation	
Annotations Comparison Quantification • Options Track water • Add track • Options Track nume • Add track • Spot information Track up §pot information Track up • Spot information • Spot information • Spot information • Spot information • Options	Plate image Densitogram	+	Plate image Densi	togram Annota	ations Comparison Quantifica	ation
Color Plate name Evaluation Color Plate name Evaluat		Annotations Comparison Quantification		~ 01	ptions	
+ - Color Plate name Evaluation + - Color Plate name - Co		Voptions Trackwidtb(mm)	REAREARE		[mm] 8	+
Color Track name thannel Color Track SWhite Original: ⊖ Spot Information Track ID		+			Color Plate name Track name Evaluation chann	nel
∧ Spot Information DyeMix 20x10 cm: Track 5 White Original: ⊖ Select a track. Select a track. Track ID Spot Information ~ Select a track.		Color Track name channel Add track			← Cm Track 5 Original ~	Ô
Spot information → Spot information → and channel Select a channel		 Spot information 	DyeMix 20x10 cm : Track 5 White Original : ⊖	Select a	Select a track 🗸	
Spot text hRf		Track ID Spot information ~ Spot text hRf		and ch	hannel Select a channel	
+ Add spot		+ Add spot			Done	
Capture/Load Image Plate Export Report		Capture / Load Image Plate Export Report		~ Sp	oot information	

图4.15:为绘制密度图选择一个轨迹







图4.17:为密度图和板上的斑点命名

3

4

5

6

7

8

9

10

11

- 在创建评估之前,已在板页面中定义了所有相关的轨迹(参见第4.7节)
- 您至少拥有用于测量的一个参考轨迹

您位于Evaluation(评估)页面

为估算物质的浓度,构建轨迹的斑点需与一或多个参考轨迹(即,物质的浓度已知)的相应斑点进行比较。一般来说, 比较是基于密度图中斑点峰的高度和面积进行的(第4.10.1节提供有关如何计算密度图的详细信息)。多个参考使您 能够通过回归方法更准确地进行估算。

4.11.1 准备定量评估

4.11 定量分析

前提条件:

定量分析由选择斑点、将某些斑点定义为参考并优化密度图计算构成,从而提供良好的回归基础。以下的逐步操作 示例图解了一个可行的方法,能够高效利用软件执行此任务(您当然可以找到更适合您条件的不同策略)。

- 1) 从Create new Evaluation (创建新评估) 按钮下方的下拉菜单中选择一个板。
- 2) 从板选择下方毗邻的Illumination(光照)下拉菜单选择板图像。
- 3) 因为只需使用红光、绿光或蓝光像素值,所以请从Channel (频道) 下拉菜单中选择它们。
- 4) 点击Quantification (定量) 标签页。
- 5) 如有需要,使用滑块调节轨迹宽度。
- 6) 通过点击+ Add Peak (+ 添加峰值) 来激活斑点选择。
- 7) 在预览图像中选择一个要进行分析和放大的斑点,方法是在触控屏上使用张合手势或在台式机电脑上使用鼠 标滚轮来操作斑点。
- 8) 点击此斑点的中心;它将被添加到Quantification(定量)标签页下方的表格中。
- 9) 视需要在表格中调节前沿峰和后沿峰的R_f值。如果您没有看到front(前沿)和end(后沿)栏位,请打开相应轨迹 的下拉菜单。
- 10) 缩小,并点击您想要包含在分析中的所有其他轨迹。软件将识别轨迹,并自动将后续其他相同的峰值位置认定 为与您添加的首个此峰值位置相同。
- 11) 缩放每个斑点(参见7)并视需要在表格中调节其前沿峰和后沿峰的R_F值。
- 12) 如果您尚未在Plate(板)页面中这样做(参见第4.7节),请立即定义哪些轨迹作为参考。
- 13) 接下来,为每个轨迹输入相应的浓度和样品数量。
- 14) 点击预览图像上方的Densitogram (密度图) 标签页。

- 15) 在斑点上方滑动手指(或按下鼠标左键)放大所选斑点,在左右两边各留下峰宽度三分之一的空白。
- 16) 展开Quantification (定量)标签页下方的Options (选项)区。
- 17) 为Densitogram Calculation (密度图计算)选择Average (平均值)或Median (中位数)。
- 18) 勾选Subtract Baseline (减去基线)复选框以便自动检测和减去基线,即,没有斑点的板的背景级别 (通过每个 轨迹左侧和右侧的空白区域在本地进行估算)。
- 19) 点击 'Savitzky-Golay' 方框执行平滑功能,并选择适合的过滤器筛选范围。
- 20) 回看表格,您会看到Offset(补偿)列(如果没有,请从表格标题下拉菜单中选择Track Data(轨迹数据))。通过更改补偿值,您可以上下移动每个峰,直至轨迹基线基本上为零。在密度图中峰以外的区域点击,会显示标尺和该位置上每个轨迹的高度-这可以较好地估算最佳补偿值。
- 21) 尽管这对实际计算没有影响,但您可能需要在轨迹输入表格上方的文本栏位中输入正确的浓度单位。这个单位 后期将显示在您能为此Evaluation(评估)而生成的导出和报告中。

要在这个密度图中看到一个或多个选定的轨迹,请展开表格中的轨迹选项(如果未展开轨迹选项,软件将绘制所有 轨迹)。在对所有峰的密度图感到满意后,您可以继续评估峰值测量和优化回归设置。

2

1

4

9

10

11



4.11.2 读取定量测量结果和微调

在板预览的下方(或,根据所选标签页,也可能是密度图下方),您可以找到总结了分析结果的两个表格,每当修改了 输入值,分析结果都会自动更新。请注意,只有为所选回归方法,正确准备了密度图和充足数量的参考轨迹,才会获 得正确的分析结果(参见前文章节)。



图4.19:选择用于定量分析的回归方法

上方表格列出了根据峰高度和面积计算得到的所有非参考轨迹的浓度。此表格末尾给出了准确的回归公式及回归 系数 'r'和标准偏差 'CV'。选择预览图像 (或密度图) 上方的Calibration Curve (校准曲线) 标签页,将显示相应的图 形,有助您评估分析的质量。根据可用参考轨迹的数量,您可以测试不同的回归方法,从中找出能够给出最准确结果 的方法。从Quantification (定量) 标签页下Options (选项) 区的Regression (回归) 下拉菜单,能够选择可选的回归 方法 (其中,c代表浓度,x代表峰面积或高度):

Linear-1(线性-1),需要至少1个参考轨迹

c = ax

Linear-2(线性-2),需要至少2个参考轨迹

c = ax + b

Polynomial(多项式),需要至少3个参考轨迹

 $c = ax^2 + bx + c$

Michaelis Menten(米氏方程),建议至少使用3个参考轨迹

$$c = \frac{ax}{(b+x)}$$

3

5

9

10

11

12

下方表格概括了用于计算的峰详细信息。在使用上方表格提供的定量结果之前,请检查下方表格提供的数值是否可用。

如果发现不合理或不需要的值,您总是能返回到上文介绍的准备步骤,进行纠正或微调设置。只需点击几下,您甚 至能够检验不同的板图像(Illumination(光照))或频道是否可以提高分析质量。两个结果表格将会因应所做变更 而自动更新,给您即时反馈。

4.12 导出图像和评估结果

前提条件:

- · 您位于Plate(板)或Evaluation(评估)页面
- 如果在相应表格中输入数据,则仅会生成某些特定的导出。

TLC Explorer本身不应是板记录的最终目的地。特别是,长期存储您的结果并不是它的设计用途。在很多应用案例中,您可能只想执行上述工作流程,直至产生需要的结果,并从设备导出结果和加以使用,这可能只是一个简单的板图像,也可能是含有密度图数据的表格,或是定量分析的详细结果。

例如,如果只想获得紫外光照射下的TLC色谱板的图像,那么您只需要按下Plate(板)页面中的Download image (下载图像)按钮。板页面会在捕捉了所有图像后自动打开。这个按钮是保存预览区中图像的快捷方式。要保存其他 图像,须通过Plate(板)、Illumination(光照)和Channel(频道)下拉菜单来配置预览,或是点击Show Overview(显 示一览)获得整个底板的图像。



图4.20:评估的导出注释对话框

7

9

11

由于Evaluation (评估)页面上的工具可生成多种数据 – 如密度图、绘制和表格 – 该页面底部的Export (导出) 按钮 会打开一个对话框,让您能够选择要导出的特定数据 (参见图4.20)。显示的对话框根据您打开的标签页而有所不同:

4.12.1 '导出原始裁切图像'(所有标签页)

这个导出包,包括在Plate(板)页面生成的数据:

- 经过裁切且底部边缘已水平对齐的所有板图像
- · 每个板都具有一个名为 'Plate_<plate name>.csv' 的文件, 它用Info (信息) 区下添加的元数据描述了板, 文件中 包括的表格则提供了所有板相关图像文件的详细信息, 例如, 捕捉图像的用户、时间和日期及曝光时间。
- · 文件 'Plate_<plate name>_Tracks.csv',包含Plate(板)页面中轨迹表格的内容。

4.12.2 '导出显示的图像'(所有标签页)

选择此导出包,能获得带有当前所选图层(即您在预览区看到的图像)的所有板图像,有关详细信息,请参见第4.8 节。唯一的例外:注释和斑点索引编号将被其相应的注释文本或斑点名称所取代。为区分这些图像和原始板图像,这 些图像的名称带有 '_overlay.jpg' 后缀。

4.12.3 '导出注释'(仅'注释')

每个板的Annotation(注释)标签页下的上方表格(参见第4.9节),会被导出到名称带有'_Annotations.csv'后缀的 文件中。除注释文本外,该文件还包括距离板左下角有几毫米远的准确位置,以及图像像素坐标。请注意,图像像素 的索引通常像一个矩阵,也就是说,原点位于左上角,后面的像素坐标直接在横纵轴上做加法。

4

11

4.12.4 '导出斑点'('注释' 和'比较')

在Comparison (比较)标签页的下方表格中命名的斑点被按照注释导出到带 '_Spots.csv' 后缀的文件中。准确位置的给出方式有两种,一种是距离板左下角几毫米远,另一种是图像像素坐标。除注释表格外,斑点包括有关其所属轨迹的信息。

4.12.5 '导出比较'('比较'和'定量')

每个用于比较的密度图,或者说是作为定量分析的基础,都可以导出为一份单独的CSV文件。出于可追溯性的考虑, 它始终包括标题为RAW_data(原始_数据)的列,此列包括了每个元素的未经平滑或基线减法的原始数据;第二列, 则是经过处理的轨迹。

要简化图形绘制或导出数据的处理,此表格中的x坐标将以三种形式提供:像素、hR_F和距离板底部的距离(毫米)。

4.12.6 '导出定量'(仅'定量')

选择此导出包,将包括定量结果文件,以及包含底层数据和所选计算设置的文件。为使屡次分析都能获得相同结果,导出应包括密度图数据和使用的板图像。

4.12.7 下载或保存导出的文件

导出对话框提供用于获取文件的两个选项。您可以把已经打包和压缩为zip归档文件的所选文件,下载到您用来控制用户界面的设备中,或将相同的zip文件保存到插接到设备背面某个端口上的USB存储棒中(B2或B3)。

2

3

4

5

6

8

9

4.13 生成有关评估的报告

前提条件:

- · 您位于Evaluation(评估)页面
- ·报告基于特定的评估。在按下Report(报告)按钮之前,请检查从区域顶部下拉菜单选择了正确的评估。

点击Evaluation (评估)页面底部的Report (报告) 按钮,让软件编译出概括分析。根据默认设置,它包含您输入的所有数据以及根据您的输入生成的所有结果。通过取消选择您想排除的区域右侧的复选框,可减少分析包含的内容。



图4.21:创建和配置有关评估的报告

展开不同区域,允许您添加诸如报告标题等更多信息,或微调您想包括在报告中的内容选择。某些栏位(如,图像名称),纯粹只为提供信息,但出于维护分析可追溯性的缘故,不能更改。

每份报告还包括序号等用于识别设备的信息和软件版本。为在日后确认在创建报告后,没有图像文件或分析数据被 更改,报告列出了这些文件的哈希值(利用MD5信息摘要算法而生成)。这类似于指纹加密。对这些文件的任何更改, 都会导致值的变化。

点击Print(打印)按钮即可将报告导出为PDF文件,然后保存下来供日后参考。请注意,PDF中的图像经过缩放,分辨率比原始文件要低。而使用File Manager(文件管理器)(参见第4.14节)或Export(导出)功能(参见第4.12.1节和第4.12.2节),可获得较高分辨率的图像。



10

12

版本1.0-2024/02

3

4

5

6

7

8

9

10

11

12

M TLC EXPLORER 10:04 PM 🗮 **Files** Only my files Search... Export Сору Delete Import Image Set Username Date ~ ~ DyeMixes EgonStahl 21/10/2023, 08:17:44 White UV366 UV254 Evaluation_1 Evaluation 2 DyeMix lot 42 EgonStahl 21/10/2023.08:34:32 < □ Remaining free space: 84.1% / 17.6GB Back 图4.22:文件管理器

4.14 管理文件和图像集

前提条件:

- · 您或其他用户已创建至少一个图像集
- · 点击主菜单图标(位于右上角的三个横条)打开文件管理器,然后选择File Manager(文件管理器)。

通过右上角的主菜单,可在工作流程的任何时点,转至文件管理器。如果要返回到从其打开文件管理器的页面,请按下页面底部的返回按钮。

所有文件和文件管理器,都通过图像集关联在一起。属于某位用户的多个图像集,包含了该用户捕捉的不同光照下的所有图像,及由这些图像衍生而得的所有板图像和数据。一个图像集可以包含多个评估,每个评估都具有作为评估依据(如,轨迹比较和定量分析)的单独一份板图像和轨迹数据,有关详细信息,请参见第4.8节。

在打开文件管理器后,表格将列出属于当前登录用户的图像集。要查看其他用户的图像集,须取消选择右上角的 Only my files (仅我的文件)复选框。该框左侧是Search (搜索)文本输入栏位,使您能够筛选出包括您输入文本的图 像集名称列表。

在文件管理器屏幕右下角,应显示Remaining free space (剩余空闲空间)。请至少保证250 MB的空闲空间,以确保 系统平稳运行。

4.14.1 导出图像集和下载图像

一项通用任务是下载特定的板图像。展开表格中的相应图像集(点击行首的箭头),是下载的快捷方式。为获得相机 拍摄的原始图像,请点击所需的显示为绿色的光照类型,此时会显示预览,您可以下载图像了。展开Evaluation(评 估)并选择光照类型,就能找到经裁切和对齐后的板图像。如果在多个原始图像中找到多个板,板图像将列在对应的 板名称下。

通常导出一个或多个选定的图像集会有多个目的,这是因为生成的zip文件包括在TLC Explorer中存储的所有相关 文件。关于CSV文件中的表格数据等结构化数据,请使用Evaluation(评估)页面中的导出选项(参见第4.12节)或是 生成一份报告(参见第4.13节)。导出的图像集主要用于在合适的外部位置上长期存储数据,同样,日后它们可以被 (重新)导入到同一台TLC Explorer设备,或传输到其他TLC Explorer设备上。

要导出图像集,须勾选所需图像集左侧的复选框,然后点击Export(导出)。打开的对话框提供了一个选项,方便您在成功导出数据后,删除存储在TLC Explorer上的数据。您还可决定将数据保存在USB存储棒上(连接到B2或B3),而 非根据默认设置下载zip文件。建议检查导出的数据,确认选择了正确的图像集。为检测可能出问题的导出,只需把 导出的文件导入(见下文,无需实际导入它):导入功能将自动检查图像集是否完整和有效。

Export Image Sets Image Set Username DyeMixes EgonStahl DyeMix lot 42 EgonStahl DyeMix lot 42 EgonStahl Export Image Sets to USB stick No USB stick found Delete Image Sets after successful export	Export Image Set Username DyeMixes EgonStahl DyeMix lot 42 Export Image Sets to USB stick No USB stick found Delete Image Sets after successful export Cancel	Export Image Set Username DyeMixes EgonStahl DyeMix lot 42 Export Image Sets to USB stick No USB stick found Delete Image Sets after successful export Cancel Export	Image Set Username DyeMixes EgonStahl DyeMix lot 42 EgonStahl Export Image Sets to USB stick No USB stick found Delete Image Sets after successful export	
Image Set Username DyeMixes EgonStahl DyeMix lot 42 EgonStahl Export Image Sets to USB stick No USB stick found Delete Image Sets after successful export	Image SetUsernameDyeMixesEgonStahlDyeMix lot 42EgonStahlExport Image Sets to USB stickNo USB stick foundDelete Image Sets after successful exportCancelExport	Image SetUsernameDyeMixesEgonStahlDyeMix lot 42EgonStahlExport Image Sets to USB stickNo USB stick foundDelete Image Sets after successful exportCancelExport	Image SetUsernameDyeMixesEgonStahlDyeMix lot 42EgonStahlExport Image Sets to USB stickNo USB stick foundDelete Image Sets after successful export	
DyeMixes EgonStahl DyeMix lot 42 EgonStahl Export Image Sets to USB stick No USB stick found Delete Image Sets after successful export	DyeMixes EgonStahl DyeMix lot 42 EgonStahl Export Image Sets to USB stick No USB stick found Delete Image Sets after successful export Cancel	DyeMixes EgonStahl DyeMix lot 42 EgonStahl Export Image Sets to USB stick No USB stick found Delete Image Sets after successful export Cancel Export	DyeMixes EgonStahl DyeMix lot 42 EgonStahl Export Image Sets to USB stick No USB stick found Delete Image Sets after successful export	
DyeMix lot 42 EgonStahl Export Image Sets to USB stick No USB stick found Delete Image Sets after successful export Cancel Export	DyeMix lot 42 EgonStahl Export Image Sets to USB stick No USB stick found Delete Image Sets after successful export Cancel Export	DyeMix lot 42 EgonStahl Export Image Sets to USB stick No USB stick found Delete Image Sets after successful export Cancel Export	DyeMix lot 42 EgonStahl Export Image Sets to USB stick No USB stick found Delete Image Sets after successful export	
Export Image Sets to USB stick No USB stick found Delete Image Sets after successful export Cancel Export	 Export Image Sets to USB stick No USB stick found Delete Image Sets after successful export Cancel 	 Export Image Sets to USB stick No USB stick found Delete Image Sets after successful export Cancel 	Export Image Sets to USB stick No USB stick found Delete Image Sets after successful export	
No USB stick found Delete Image Sets after successful export Cancel Export	No USB stick found Delete Image Sets after successful export Cancel Export	No USB stick found Delete Image Sets after successful export Cancel Export	No USB stick found Delete Image Sets after successful export	
Delete Image Sets after successful export Cancel Export	Delete Image Sets after successful export Cancel Export	Delete Image Sets after successful export Cancel Export	Delete Image Sets after successful export	
Cancel	Cancel Export	Cancel		
			Cancel	Export

图4.23:文件管理器中的导出图像集对话框

1

3

6

12

4.14.2 复制或导入图像集

用户只能加载和处理他们自己的图像集。要操作其他用户的图像集,请取消选择右上角的Only my files (仅我的文件)选项,查找和选择所需的图像集,然后点击Copy (复制)来创建属于当前登录用户的一份拷贝。您当然可以复制您自己的图像集。

也可在同一台或其他多台设备上导入被导出的图像集,这不仅允许恢复数据,而且还支持您与其他使用 TLC Explorer的实验室共享图像集。数据须为原始zip格式、完整和无损坏,否则导入功能将拒绝导入它。

		×	
e Sets			
Overwrite?	Import as		
	Herbal extracts		
ing Image Sets			
Import			
	CVERWRITE?	CVerwrite? Import as Herbal extracts ing Image Sets Import	e Sets Vverwrite? Import as Herbal extracts ing Image Sets Import

图4.24:文件管理器中的导入图像集对话框

4.14.3 删除图像集

建议您定期检查在文件管理器右下角显示的剩余存储容量,并删除过时的图像集。

在删除数据之前,您可以备份整个设备(参见第6.1节),或将所选图像集导出到合适的长期存储位置(参见 第4.14.1节)。

当您确定不再需要存储在TLC Explorer上的图像集后(例如,已在别处存储了拷贝),可勾选左侧的复选框,然后 点击Delete (删除)。系统将在永久删除数据之前,要求您确认确实删除列出的图像集。此操作无法撤消。

4.15 重置或关闭设备

和所有台式机电脑一样,在拔下电源插头之前,需要正确关闭TLC Explorer。否则,可能损坏设备。 您可以按照第3.6节"关闭设备"中的介绍,查找通过不同方式和选项来重启或关闭设备的详细信息。

11

5 用户管理和设备设置

前提条件:

· 您位于Settings(设置)页面。通过主菜单中的Settings(设置)选项,可抵达该页(参见图4.3)。

	TLC EXPLORER	10
Management	Network	Illumination settings
User management	Device name	Stray light detection
Backun data	tlc-explorer-da230010	Apply illumination correction
Duckap data	Device serial number	Save
Restore backup	da230010	Create illumination correction data
	Pacaive IP from DHCP for Ethernat	
	Current Ethernet IP	
Update from USB drive	192.192.192.210/24	
Recover old version		
Factory reset	Provide own WiFi network	
Current version : 01.01.40-RC10	tlc-explorer-da230010	
Report template	Password	
generator		
Upload brand image	192.168.200.1/24	
Upload CSS format template		
	Save	

图5.1:设置页面

- 多位用户可同时在TLC Explorer上工作。这使不同用户能够访问和处理此前记录的板图像,与此同时,另一位用户则可捕捉新的板图像。然而,实际上一次只能获取一个图像集中的多幅图像,系统会协调多个并发活动,或者说是 "会话",以确保满足下列两个条件:
- · 仅一位用户可以捕捉图像。
- 一位用户仅可在一个浏览器中登录(例如,在平板电脑或笔记本电脑上)。
- 如果第二位用户尝试在其他用户正在捕捉图像的时候也去捕捉图像,系统将发出警告并锁定此功能,直至第一位用 户完成了图像捕捉过程。
 - 用户在欢迎屏幕上登录(参见图3.5),将自动终结该用户在其他多台设备上的所有其他会话。

2

3

4

5

6

8

9

10

11

5.1 更改活动的用户

打开主菜单后,选择Change User(更改用户)选项,即可更改活动的用户(如图4.3所示)。这将注销您的当前用户身份,将您重定向至登录屏幕(图3.5),在这里可以选择新用户,然后按下Login(登录)按钮。

5.2 添加和删除用户

从主菜单选择Settings(设置),然后点击User Management(用户管理)按钮,就可以添加或删除用户。 要添加新用户,请点击Add a new user(添加新用户),然后在显示的对话框中输入用户名和全名(全名是可选项,例 如,在报告中使用)及用户语言偏好。

要删除用户,可点击列出所有用户的表格中的回收站图标即可。屏幕上会打开一个对话框,要求确认是否删除,并且还会显示一个复选框,供您选择是否同时删除与用户名相关的数据。如果在点击Delete (删除)之前未勾选这个复选框,则该用户的图像集将保留在设备上,可在文件管理器中用户名 'DELETED_USERS' 下找到 (参见第4.14节)。



图5.2:添加和删除用户

11

1

2

3

4

5

6

7

8

9

5.3 设备名和序号

在Network (网络) 标题下, 您可以找到Device name (设备名) 和Serial number (序号)。后者不能更改, 且应与您在 设备外壳上看到的数值一致 (见B6)。

用户可以设置Device name(设备名),设备名将用于在报告、导出的数据和网络中识别设备(参见第8.2节)。尤其是 在您实验室部署了多台TLC Explorer的情况下,独一无二的名称将帮助您由结果(如,导出的板图像)逆向追溯回至 正确的设备。请牢记,按下Save(保存)按钮才能应用您做的更改。

5.4 时间和日期

设备关闭后,TLC Explorer没有内部时钟运行。因此,软件会尝试从笔记本电脑或平板电脑等用户设备获取当前的时间和日期。另一个可选方法是在Setting(设置)页面手动设置日期和时间。请牢记,按下Save(保存)按钮才能应用您做的更改。

5.5 数字光照校正

要改进光照的均匀性,您可以根据来自参考板的一系列图像,在电子文档上对图像进行数字化校正。 要校准系统,请将一张白色、平整、20.5厘米 x 20.5厘米大小的复印/打印纸放置在底板上并关闭抽屉。然后,按下 Illumination Settings(光照设置)标题下的Create illumination correction data (创建光照校正数据)按钮,并按 照屏幕上的指示进行操作。切勿在校正过程完成之前,打开抽屉或检测窗。系统会在所有可用光源的照射下拍摄多 幅图像,然后为每种光照类型计算校正表。

请注意,错误校准的光照校正可能导致图像带有伪影,降低结果的质量。要检查校准是否带来了希望的改进,请将底 板上的参考板手动旋转90度,捕捉测试图像集。忽略每个边缘处约1厘米的空白,从各个角度观查板应具有同等亮 度。

您可以通过Apply illumination correction (应用光照校正)复选框来激活或停用数字光照校正。

5.6 杂散光检测

激活Stray-light detection (杂散光检测)后,系统会在捕捉图像之前检查检测窗是否正确关闭。如果在无光源开启的情况下底板不是全黑,则软件会发出错误信息并停止捕捉。

建议您启用此功能,它有助于防止捕捉的图像带伪影或是对比度下降,因此提高了所有后续处理步骤的数据质量。

4

6

9

10

11

系统提供多个选项来导出在TLC Explorer上生成的数据。如果您希望保存精选的图像集,使用文件管理器更快速更 灵活(参见第4.14节)。备份功能将获取设备上的所有用户数据(不仅是当前登录用户的数据),并将其压缩为一个文件,它可用于在同一台设备上恢复用户设置或是在其他TLC Explorer上克隆这些用户数据。

6.1 生成备份

从主菜单选择Settings(设置),然后点击Management(管理)标题下的Backupdata(备份数据)。在显示的对话框中(参见图6.1),可选择是通过网页浏览器下载备份文件,还是将备份文件存储在连接的USB棒上。

根据默认设置,备份包括所有图像集,但不包括用户设置。为了在备份中包括用户设置,请在选择存储选项之前勾选 相应的复选框。

备份可能需要数分钟才能完成,这取决于您在设备上存储的图像集和评估的数量。

 ×
Backup data
Include settings: users, device settings, illumination correction data
Download with browser Backup to USB

图6.1:备份设备上所有用户数据和设置的对话框

10

11

6.2 恢复备份

同一个对话框(在点击Management(管理)标题下的Backup data(备份数据)后打开),提供通过USB驱动器或是 上传备份文件的方式来恢复备份。一旦用户确认后,备份数据将覆写所有用户数据和设置,系统可能需要数分钟来 完成这个过程。然后系统将自动重启,当设备准备好再次操作后,浏览器应导航至登录页面。

6.3 将用户数据迁移至不同的TLC Explorer

在不同设备上克隆用户数据和设置的方法,与在同一台设备上恢复备份数据的方法相同。但是,由于目标设备上不同的硬件和可能不同的软件版本,系统会提示两条警告信息:

- · 要确保图像集数据的兼容性,请确保目标设备运行与原始备份设备相同或更新的软件版本。理想做法是在开始 恢复过程之前,更新至最新软件版本。
- 备份可能包括光照校正数据。很可能出现的情景是,按照第5.5节描述的步骤为新设备重新计算校正表。

2

3

4

5

6

8

9

10

11

7 软件更新

前提条件:

· 您位于Settings(设置)页面。通过主菜单中的Settings(设置)选项,可抵达该页(参见图4.3)。



7.1 当前软件版本

您可在Management (管理) 标题下的Settings (设置) 页面中看到当前软件版本,最后一行显示了准确的版本号 (参见图7.1)。

7.2 获取软件更新

为使TLC Explorer使用新功能和安全性改进而保持最新状态,请在TLC服务网站上定期检查更新的软件版本。通过 扫描图7.2中显示的二维码或在网页浏览器中手动输入链接https://sigmaaldrich.com/TLCservice,您将被定向至 提供TLC Explorer产品和可用软件更新的最新信息的网站。

只需将该网站上的最新软件更新文件下载到您的电脑或移动设备上即可。

产品ID 1.52610.0001明确标识出了兼容的软件。

请注意,更新须由用户发起,即使在TLC Explorer接入互联网时,它也不会自动搜索更新。



图7.2:软件更新位于:https://sigmaaldrich.com/TLCservice

7.3 安装更新

系统提供将软件更新传输至您的TLC Explorer的两个选项:您可以在网页浏览器中直接上传文件,或将该存储棒连 接至设备然后复制USB存储棒上的更新文件。

通过网页浏览器上传更新是更简便和推荐的方法:在Management (管理)标题下的设置页面中,点击 Upload update (上传更新),并从您的电脑或移动设备中选择文件,然后点击Upload update (上传更新)按钮。 更新过程会耗时数分钟,其间包括重启TLC Explorer。切勿在更新时断开设备电源并等待更新过程结束。一般来说, 浏览器将自动识别TLC Explorer是否可以重新操作,然后将您重定向至Login (登录)页面。如果浏览器不支持这一 机制,只需花费一分钟等待电源图标F1停止闪烁并以明亮白光亮起后,打开TLC Explorer网站。

要从USB存储棒更新TLC Explorer,请复制该存储棒根目录中的更新文件,也就是说,更新文件不位于文件夹中。当 将存储棒连接至USB-A端口B2或B3后,请点击Update from USB drive(从USB驱动器更新)(参见图7.1)。屏幕上 会显示一个对话框,让您选择想要安装的软件版本(此时,存储棒上应有多个更新文件)。请让系统花费几分钟完成 更新过程。

如果更新过程发生错误,系统将自动回退至启动更新时所安装的软件版本。

3

6

11

4

5

6

7

8

9

10

11

12

在TLC Explorer的启动过程中(即,在电源图标F1闪烁的过程中),触控并按住全部三个光照按钮

置网络设置,请参考第8.4节。

(F2、F3和F4)。

7.5

1.

2.

7.4 恢复前一软件版本 用户可视需要手动回退至安装的前·

恢复出厂设置

系统提供两个启动恢复出厂设置的选项:

用户可视需要手动回退至安装的前一软件版本。此时,点击Recover old version (恢复旧版本)按钮,会打开软件降级菜单并列出可用的旧软件版本。系统只会存储一个旧版本以及设备交付时安装的软件版本。只需选择所需版本,然后点击按钮安装它。与更新过程相同,降级过程会耗时数分钟,其间包括重启TLC Explorer。如果您希望安装的旧版本与软件降级菜单列出的版本不相同,您可以使用前文介绍的安装过程(参见第7.3节)。

发起恢复出厂设置应格外谨慎小心,这会擦除所有用户及其数据,以及包括网络设置在内的所有设置。如果只需重

点击设置页面中Management (管理)下的Factory reset (恢复出厂设置),在打开的对话框中确认您的意图。

恢复出厂设置还会重新安装设备交付时所安装的软件版本,而擦除设备中所有其他软件版本。

8 连接TLC Explorer

TLC Explorer具备多种连接选项,可轻松集成到实验室的现有网络中,且提供它自己的联网基础设施。更多设置选项见下文。然而,将设备集成到您的实验室网络中,可能会有意外后果。根据网络配置,访问网络的每个人都可以访问TLC Explorer设备。用户数据或设备设置都不受密码保护,这是因为假定实际的或是数字访问都由实验室经理控制。

请注意:仅应在咨询您的网络管理员后,才能将设备连接到现有网络中。

初次连接您的TLC Explorer,请参考第3.4节,其中对把笔记本电脑或平板电脑链接至TLC Explorer和打开用户界面 有详细介绍。

Device name	
tic-explorer-da230010	
Device serial number	
da230010	
Receive IP from DHCP for Ethernet	
Current Ethernet IP	
192.192.192.210/24	
Provide own WiFi network Network name (SSID) tlc-explorer-da230010	
Password	
Current WiFi IP 192.168.200.1/24	
Savo	

图8.1:在设置页面的"Network(网络)"区配置连接设置。

8.1 TLC Explorer自己的WiFi

只要没有更改出厂设置或在恢复出厂设置(参见第8.4和7.5节)后,设备会广播它自己的WiFi信号。通过扫描出厂交付包装中的贴纸上的二维码,可以轻松连接到此网络中。二维码1包含平板电脑或智能手机连接至TLC Explorer自有WiFi网络所需要的全部信息。扫描并按屏幕上显示的指示进行操作。

另一种可选方式是,可以在平板电脑或具备WLAN适配器的台式机电脑上打开网络设置,搜索以'tlc-explorer'开头 且后跟序号B6的网络名称(SSID)。选择'WPA/WPA2'为安全性设置,输入二维码贴纸上的默认密码("步骤1"中设备 名后面就是密码)建立连接。

移动设备可能警告您所选WiFi没有互联网连接,根据设计这属于正常情况。但是,某些移动设备会首选具有互联网可访问性的网络连接。因此,有必要打开移动设备的WiFi设置并禁用具有互联网可访问性的已保存网络的Auto-connect(自动连接)-除TLC Explorer外,这些已保存网络的互联网连接也是可用的。

如果平板电脑、笔记本电脑或智能手机是首选用于控制TLC Explorer的设备,我们建议激活TLC Explorer WiFi的自动连接,并为所有其他网络停用此WiFi连接。否则,您将需要在每次网络连接中断后,手动重新选择TLC Explorer的这一网络(如,因为重启平板电脑或TLC Explorer后)。

您可以在设置页面的Network(网络)区下,更改TLC Explorer广播的WiFi网络名称(SSID)。在那里,还可以设置新的WiFi密码,用以保护网络连接和设备。在按下Save(保存)以确认更改之前,请确认勾选了Provide own WiFinetwork(提供自己的WiFi网络)前面的复选框。

8.2 通过以太网线缆集成到LAN中

在将以太网线缆插接到RJ45端口B4之前,请先联系网络管理员。根据默认设置,将从网络中的DHCP服务器为 TLC Explorer分配一个IP地址。如有必要使用静态IP地址,请取消选择Receive IP from DHCP for Ethernet (接受 DHCP分配的以太网连接IP地址)前面的复选框,并在Current Ethernet IP (当前以太网IP地址)栏位中输入静态IP 地址。要再次接受DHCP服务器分配的IP地址,将需要勾选此复选框,点击Save (保存)然后重启设备。 如果连接的LAN具有本地DNS服务器,您可以从网络中的任何其他设备连接至TLC Explorer,方法是在设置页面中 的Network (网络)下,输入后接设备名的http://。尤其是如果您实验室使用多台TLC Explorer,请为其分配独一无 二的设备名,然后在网页浏览器中把基于地址的相应设备名加为书签 (例如,http://tlc-explorer-quality-lab,或 http://tlc-room214)。如果没有可用的本地DNS,还可以通过Current Ethernet IP (当前以太网IP地址)标签页下显 示的IP地址访问每台TLC Explorer,比如,在浏览器地址栏中输入http://192.168.178.210 (无'24')。 3

2

4

9

8.3 集成到现有WiFi网络中

TLC Explorer还可以连接到现有WiFi网络中,而非提供自己的WiFi网络。然而,如果可以通过以太网线缆连接设备,则此选项通常根据用户偏好设置。

要设置与现有WiFi网络的连接,需要准确的网络名称(SSID)和该网络的密码。取消选择Provide own WiFi network (提供自己的WiFi网络)前面的复选框,在下方的栏位中输入网络名称和密码。仔细检查是否正确输入了名 称和密码,输入错误会导致无法访问TLC Explorer,并需要按照第8.4节的描述重置网络连接。只有确认输入正确后, 才能点击Save (保存)。

与通过以太网连接类似(参见第8.2节),只需在网页浏览器地址栏中输入后接设备名的http://(如果有可用的DNS) 或设备IP,便可从WiFi网络中的任何设备访问TLC Explorer。

8.4 重置网络连接

如果您所做的网络设备更改,或是遗忘了WiFi密码,妨碍您连接至TLC Explorer,您总是可以将网络设置恢复为出 厂默认设置。要重置网络连接,您需要按下电源图标F1直至图标背光变为暗淡白光,关闭设备。在TLC Explorer重启 过程中,触控并按住全部两个光照按钮(F3和F4)。重启过程是指电源图标F1闪烁的过程。切勿按下全部三个光照按 钮。这将引发彻底恢复出厂设置,删除所有用户数据和设置(参见第7.5节)。

8.5 将TLC Explorer作为网络位置(SMB)安装

除通过文件管理器导出图像和图像集外(参见第4.14.1节),也可以通过运行在TLC Explorer上的samba服务器访问 所有用户数据。与TLC Explorer在同一网络中的Windows用户,应可在Windows资源管理器中Networks(网络)下 找到列出的设备名,而Mac用户应在Finder(查找器)窗口下看到设备名。Ubuntu用户需要在Nautilus窗口中按下 Ctrl+ 'l',然后输入后接设备名的smb://(如果有可用的DNS)或设备IP地址。不受操作系统的影响,访问不要求任 何用户名或密码(匿名),并应显示名为data(数据)的文件夹。此文件夹下包含为每个用户建立的属于他的图像集子 文件夹。

为保护数据完好性,用户可以读取和复制您本地台式机或笔记本电脑上的图像集,但不可以在此图像集目录上写 入。

3

4

6

8

11

10

9 故障排除

识别错误 9.1

在前面操作中您未发现的错误,有时会导致系统提示意外行为或错误。要加速搜索以查找错误的出处和潜在解决方 案,图9.1列出了一系列的检查。一旦找到问题,请执行可行的补救措施,它们显示在问题的右侧并带有编号。 切勿在拔下系统电源之前,未正确关闭设备(参见第3.6节)。这样做可能严重损伤TLC Explorer上安装的电脑系统, 造成其他问题。关闭和重启设备的安全的替代方法,是按住电源按钮F1。



图9.1:识别问题,找到解决方法

2

3

10

9.2 可行的补救措施和解决方案

- 要重启设备,用手指轻按电源图标F1,直至图标开始闪烁(指示设备正在关闭)。图标停止闪烁后维持暗淡光芒, 请等待几秒钟,然后再次按下图标重启设备。
- 2) 如果电源图标不再亮起,设备可能未连接到有效的电源。检查电源的简单方法,是将平板电脑或手机连接到设备背面USB-C端口上插接的USB-C连接器。但是,这个测试仅适用于推荐给TLC Explorer的USB-C电源适配器(参见第2.4节)-第三方电源适配器和线缆可以给手机充电,但在规格上与TLC Explorer不兼容。如果电源适配器和线缆能够为除TLC Explorer外的其他设备供电,您可能需要联系设备销售商寻求技术支持(参见第10.2节)。
- 3) 如果抽屉未完全关闭,请确认底板下方的对准销P3插入了抽屉底座P4,以便底板平稳就位于抽屉中(参见第 2.3.3节)。不应在底板上放置高于底板8 mm的物体(参见第2.3.1节)。如果抽屉似乎被设备中的物体堵住(阻 挡),您可以打开背板B7并取出障碍物。
- 4) 无论何时,只要传感器指示抽屉未正确关闭,红色的抽屉打开指示灯F6就会亮起。抽屉打开后,无法打开光源和 拍摄图像。检查抽屉是否关闭正确。偶尔,软件或传感器问题会导致系统无法正确检测抽屉是否打开。如果您怀 疑是这种情况,请重启设备。
- 5) 出于安全的考虑,在抽屉正确关闭和背板B7正确安装的情况下,TLC Explorer仅允许打开光源。同时,在另一个 用户捕捉图像时,系统会停用手动光源开关F2、F3和F4。
- 6) 如果一个或全部光源不工作,或者照射时一侧是黑暗的,则可能需要更换光照装置(参见第10.4节)。如果更换的光照装置照射时仍然是黑暗的,请确认是否插接了装置的电源连接器(参见图10.3)以及是否正确安装了背板B7。
- 7) 经由以太网连接到设备,也能解决WiFi配置问题(参见第8.2节)。可通过以太网连接访问设置页面并纠正配置 (参见第8节)。
- 8) 要将所有网络设置恢复默认值,请按照第8.4节中的说明进行操作。然后,按照第3.4节中说明的步骤进行操作, 应可连接到TLC Explorer自己的WiFi网络。
- 9) 网络管理员可能已经安装了有关网络配置或您设备设置的安全措施,这阻止了您设备连接到TLC Explorer。如 有任何疑问,请联系您的网络管理员。
- 10) 如第8.1节所述,某些移动设备会优先选择具有互联网访问能力的网络连接,TLC Explorers自己的WiFi网络 不具有互联网可访问性。因此,有必要打开移动设备的WiFi设置并禁用具有互联网可访问性的已保存网络的 Auto-connect(自动连接),这是因为除TLC Explorer外,这些已保存网络的互联网连接也是可用的。
- 11) 请注意,每台TLC Explorer广播的WiFi信号都具有不同的名称(通常包括它的序号),也很有可能带有自己的密码保护。如果您的实验室部署了多台TLC Explorer设备,则可使用各设备的贴纸上的二维码进行连接(只要默认网络设置不曾更改)。
- 12) TLC Explorer会尝试通知同一网络中的所有设备,可使用后接设备名的URL "http://"连接到 TLC Explorer(参见第5.3节)。然而,某些网络可能不支持此功能,需要用户通过它的IP地址解析设备,例 如,"http://192.168.200.1"。如果多台并发的移动设备使用同一个移动数据互联网连接,也许有必要尝试在万 维网中搜索后接设备名的URL "http://"。
- 13) 有些浏览器会强行建立加密连接,可能会自动为输入的地址加上"https://"前缀,然后会生成提示网站可能不 安全的警告信息。浏览器无法知道您正在访问受保护局域网中的TLC Explorer,其风险远比从互联网打开网站 小得多。另外一种方法是,在浏览器设置中禁用https-only(仅https),另请参见第3.4节。
- 14) Capture Now (立即捕捉)按钮仅在下列情况下启用:如果输入了独一无二且有效的图像集名称,并且系统未从 事图像捕捉。如果用户多次按下位于Use manual Exposure (使用手动曝光)区的Auto (自动)按钮来寻找最佳 曝光时间,可能会发生超长的延时。此过程可能需要一点儿时间,尤其是在使用紫外线光源照射时,这是因为必 须捕捉和处理一系列图像。

3

8

9

10

- 15)为确保没有杂散光影响图像质量,例如,由于检测窗F7打开而招致了杂散光,相机将会在开始图像捕捉之前,检查在无光源照射时底板是否全黑。建议您启用这项检查,但您仍然可以在Settings(设置)页面中禁用它(参见第5.6节)。
- 16) 同一时间仅一位用户可以捕捉图像。尽管其他用户也可加载和处理图像集,但若同时捕捉图像将会导致提示错误信息。
- 17) 如果可用存储空间不足以保存捕捉的图像,系统将会显示错误信息。请将所有当前不需要处理的数据转移至 其他设备,并在本台TLC Explorer上删除它们。在文件管理器(参见第4.14节)屏幕右下角,应显示Remaining free space(剩余空闲空间)。请至少保证250 MB的空闲空间,以确保系统平稳运行。
- 18) 软件会搜索多个这样的TLC色谱板:尺寸最小且呈明亮矩形。要使设备能够自动检测多个色谱板,放置它们时须相互间隔几毫米,以使设备能够检测到在每个色谱板周围都有至少2毫米宽的黑色涂层。
- 19) 如果您透过检测窗F7看到了正确光照下的色谱板,但是板图像上的光照不均匀,则可能已经应用了错误光照校正。您可以停用校正步骤,或重新校准计算。有关详细信息,请参考第5.5节。
- 20) 如第7.5节所述,最后的手段是通过恢复出厂设置,删除所有用户数据、设置和软件更新。
- 21) 请联系TLC Explorer的销售商寻求技术支持。

2

版本1.0-2024/02

10清洁、维护和停用

10.1 清洁

建议您使用潮湿的无绒纸巾定期清洁TLC Explorer,尤其是底板。为清除较柔软的污垢,还可以用清洁纸巾蘸取1:1 比例混合的酒精和水溶液。

在开始清洁处理之前,应先关闭设备(参见第3.6节)和断开USB-C电源B1。只有这时,才能从抽屉中取下底板,将其彻底擦拭干净。如果来自TLC色谱板的硅颗粒,或是灰尘,弄脏了抽屉或壳体,您可以使用相同的方法清洁抽屉和壳体。但请确保不要触碰光照装置。另一种可选方法是,打开背面罩板B7(参见第2.2.6节),小心地以低功率使用手持式真空吸尘器和纸巾,清洁设备内部的灰尘。

请注意:应立即清除泼溅的液体或其他物质(另请参见第1.2.1节)!

10.2 联系技术支持

有关保修、保修相关争论或设备的特定问题,请联系您的TLC Explorer销售商。同样,如果在面对第1.2.1节 '在以下情况中,切勿操作TLC Explorer...'下列出的任一情况而不确定该如何进一步处理时,请联系设备销售商或通过www.sigmaaldrich.com联系本地经销商。

有关常见技术问题和软件更新(参见第7.2节),请使用下图10.1中的链接,访问官方TLC Explorer支持网站。



图10.1:查找TLC Explorer服务网站,请至: https://sigmaaldrich.com/TLCservice

2

3

4

5

6

8

9

10

11

10.3 订购附件和更换部件

目前,TLC Explorer产品系列包括的以下产品,可作为附加设备或更换部件而单独订购:

商品编号	名称	说明
1.52610.0001	TLC Explorer	TLC记录系统的交付清单,包括第3.1节列出的部件。
1.52613.0001	光照装置	更换部件,仅当您TLC Explorer中的光源损坏后需要更换。装置内含适合所有光照类型(即,可见光以及紫外光)的LED光源和光学元件,只能整体更换。
1.52612.0001	底板	一个额外的底板可以简化和加速工作流程。当一个底板准备就绪时,另 一个底板正在TLC Explorer中进行板的记录(参见第2.3节)。
1.52611.0001	移动电源	除电网外第二种为TLC Explorer供电的方式,例如,使用设备时没有市电条件。
1.52618.0001	防尘罩	可更换的保护罩,防止TLC Explorer蒙灰。

通用TLC耗材,如TLC色谱板、溶剂和染色以及TLC处理过程所使用的工具(例如,展开室),可从Sigma Aldrich网站 www.sigmaaldrich.com订购。



图10.2:查找TLC耗材和工具,请至: https://sigmaaldrich.com

10.4 更换光照装置

TLC Explorer采用LED光源,与传统光源相比,具有超长使用寿命。特别是发射紫外光的LED光源。但它会随着时间的推移而退化,或者在某个不确定的时间点就会失效



图10.3:更换光照装置

出于安全考虑,光照装置(内含所有光照类型的LED光源)仅可整体更换。它安装在测量室上方的滑轨中。要进行更换,请按照下列步骤进行:

- 1) 如第3.6节所述关闭设备。
- 2) 转动TLC Explorer,使设备背面朝向您。
- 3) 拔下USB-C连接器B1,断开设备与电源的连接。
- 4) 用手卸下四颗螺钉,如有必要,可以使用螺丝刀取下背面罩板B7。请注意,可把螺钉套在罩板上的螺孔中,以防止丢失螺钉。
- 5) 小心拔下图10.3中标记为"电源连接器"的线缆。
- 6) 卸下光照装置左侧的固定螺钉。
- 7) 沿着滑轨拉出旧的光照装置。
- 8) 小心推动新的光照装置滑入滑轨,确保装置正确就位。请小心操作,切勿碰触玻璃组件,否则会影响光照和影像记录质量。
- 9) 重新拧紧固定螺钉,插接新装置的电源连接器。
- 10) 重新安装背面罩板,拧紧四颗螺钉。
- 切勿将旧光照装置丢弃在未分类的城市垃圾或家居废物中。相反,它应被带到电气和电子废品的官方收集处与回收 中心,或是返回至新光照装置的销售商处。

3

8

9

10.5 停用、储存和处置

如果决定停用TLC Explorer,您可以把它储存起来以便日后使用,也可以将其返回至TLC Explorer的销售商或经销商处。

TLC Explorer不具有电池。因此,您无需在储存、运输或返回设备之前,拆卸任何电池。

然而,内部存储上可能会有敏感数据。强烈建议您发起恢复出厂设置(参见第7.5节),从而在停用TLC Explorer之前, 删除所有用户数据。

还应为设备清洁掉所有可见的残留硬质物体。如果设备接触了有害化学物质或传染性/生物危害性物质,必须进行 相应地去除污染处理。请注意,如果未正确去除污染,销售商能够拒绝收集停用的设备。

要储存设备,请取出所有TLC色谱板并确认未插接USB存储棒。切勿把底板或任何附件留在设备内部,而是把底板和 附件单独打包并放置在设备旁边。理想情况是使用原始包装(包括内衬),以便在储存过程中保护设备。请确保储存 地点的干燥性,相对湿度应在20-80%之间,温度应维持在15-40°C范围内。

根据欧盟关于电气和电子废品管理(WEEE)的指引2012/19/UE,本产品使用寿命终结后,不得作为未加分类的市政废品或家居废物处置。相反,它应被带到电气和电子废品的官方收集处与回收中心,或是返回至TLC Explorer的销售商处。有关您当地回收计划的更多信息,请至https://www.sigmaaldrich.com/EN/en/life-science/quality-and-regulatory-management/recycling。





图10.4:仅把设备送至官方收集点和回收中心,或是返回给您的销售商。

11术语表

本术语表重点解释在TLC Explorer软件环境中具有特殊含义的术语。有关薄层色谱(TLC)的一般术语和解释,请参考 第1.3节。

图像集

图像集包含了用户在不同光照下捕捉的所有图像和所有板图像,以及由这些图像衍生而得的数据(参见"评估")。 对于软件而言,每个图像集都是用户数据文件系统中的一个单独的目录,在从文件管理器导出图像集后,目录的显 示将变得突出。

图像集只能由其所有者进行编辑。如果其他用户希望使用图像集数据,他们可以创建拷贝,拷贝归其所有。对复制的 图像集所做的更改,不会影响原始图像集。

评估

"评估"的英文Evaluation以大写字母E开头,是指数据结构,该结构用于存储设置和用户操作图像集而获取的结 果。用户可以从不同层面评估一个图像集里的多幅图像,并可将带有注释的每个分析、轨迹比较和/或定量分析保存 为一个单独的评估。

每个评估包括作为评估依据的板图像和轨迹数据副本。这可确保日后在板区域中所做的更改(如,轨迹数据的更改),不会影响现有评估,只会影响在更改之后所创建的评估。

板和板图像

在捕捉使用用户所选光照类型拍摄的图像后,软件会自动检测这些图像中的TLC色谱板。找到的板会被裁切、旋转并 在水平方向上对齐底部边缘,然后另存为单独的图像文件。不同光照类型拍摄的同一个板的图像,将一起存储在图 像集的一个子目录中。

在用户界面的Plate(板)页面中(参见第4.6节),您可以为每个板添加元数据,定义轨迹位置和属性。如密度图计算或 斑点测量等后续处理步骤,都基于板图像文件。

光照

Illumination (光照) 下拉菜单, 允许您选择在所需光照类型下拍摄的所选板的板图像, (例如, 可见光, 长波紫外线, 或短波紫外线, 有关详细信息, 请参见第2.1.2节)。

经过修改的板图像保存在Plate(板)页面中(例如对比度增强后的图像),会在Illumination(光照)下拉菜单中增加一个条目,该条目可用于处理原始板图像等。

轨迹基线和密度图基线

术语"基线"是指

· 起始线,将样品点样到TLC板起始线一端,或

・ 密度图的基础。

尽管轨迹基线是轨迹的(h)R_f测量的起点,绘制图形时会自动减去密度图基线,或者在设置补偿值后手动进行补偿。

2

3

4

5

6

7

8

9

10

11
12.1 设备测量和规格

(外部)电源	USB-C电源装置/ 65 W
电源	外部电源装置(HA65NM170)具有单独的电源线(1.8米长),配备了B、G、N、I、D、E/F插头 (插头必须符合本地规定)
功率要求	外部电源装置电压为100 V - 240 V,频率为50 - 60 Hz
功耗	标准工作条件:18.8W;待机模式:0.83W
测量技术	使用视频密度测量的TLC板记录系统
波长范围	TLC Explorer为检测TLC板提供标准光照(另请参见第12.2节): 可见光(VIS) 短波紫外线(UV-C) - 波长:254 nm 长波紫外线(UV-A) - 波长:366 nm
相机	RGB传感器,具有3280 × 2464有效像素; 一个像素可捕捉约85 × 85 μm的TLC板面积
通讯接口	 · USB:2个USB-A(直接连接USB棒) · 以太网:LAN · WLAN IEEE 802.11 b/g/n (2.4 GHz)
防护(IP)等级	IP2X(主机箱)
环境条件:温度	· 设备可在15°C到40°C的温度范围内工作。 · 储存和运输温度应在5°C到40°C的范围内。
环境条件:相对湿度	· 设备的工作湿度范围为20%RH到80%RH,应确保非冷凝 · 储存和运输湿度范围为15%RH到95%RH,应确保非冷凝条件
环境条件:海拔	<2000米
尺寸	374 × 312 × 290毫米(长 × 宽 × 高)
TLC色谱板尺寸	使用最大尺寸为20厘米 × 20厘米的TLC色谱板
重量	约10.4千克
仪器安全	符合 低压指令2014/35/EU IEC 61010-1, ANSI/UL EN 61010-1 CSA-C22.2 No. 61010-1-12
EMC	符合 EMC指引2014/30/EU EN/IEC 61326-1 / EN 301 489-1 FCC
RF	符合 RED 2014/53/EU EN 300 328, FCC 包含: FCC ID 2ABCB-RPICM4 IC:20953-RPICM4

12.2 光照频谱





版本1.0 - 2024/02

我们尽我们的所知与所能,向客户提供关于应用技术与法规问题的信息和建议,但恕不承担任何责任和义务。任何情况下,客户均须遵守现行法律和法规。该规定在第三方的任 何权利方面同样适用。我们提供的资料和建议并不免除我们的客户检查我们的产品是否适合预期用途的责任。

默克生命科学业务部在美国和加拿大以MilliporeSigma名称运营。

Merck Life Science KGaA, 64271 Darmstadt, Germany, +49(0)6151 72-2440

www.sigmaaldrich.com

© 2024年版权归Merck KGaA, Darmstadt, Germany和/或其附属公司所有。保留一切权利。Merck和Supelco是Merck KGaA, Darmstadt, Germany或其附属公司的商标。所有其他商标均归相应所有者所有。商标的详细信息可通过公知资源了解。

