

# Accela Open Autosampler

## Hardware Manual

60357-97102 Revision D December 2013



**Thermo**  
SCIENTIFIC

© 2013 Thermo Fisher Scientific Inc. All rights reserved.

EQuan, QuickQuan, TSQ Vantage, and Velos are trademarks, and Accela, Exactive, TSQ Quantum, and Xcalibur are registered trademarks of Thermo Fisher Scientific Inc. in the United States.

The following are registered trademarks in the United States and other countries: Agilent is a registered trademark of Agilent Technologies Inc. Cheminert, Valco, and VICI, are registered trademarks of Valco Instruments Company, Inc. Kalrez and Teflon are registered trademarks of E.I. du Pont de Nemours & Company. Microsoft, Vista, and Windows are registered trademarks of Microsoft Corporation. PAL is a trademark of CTC Analytics AG. Rheodyne and the Rheodyne logo are registered trademarks of Rheodyne, L.P. Rytan is a registered trademark of Chevron Phillips Chemical Company. Simriz is a registered trademark of Carl Freudenberg KG Corporation. Torx is a registered trademark of Textron Inc. Tygon is a registered trademark of Saint-Gobain Performance Plastics Company. Unimetrics is a registered trademark of Unimetrics Company.

PEEK is a trademark of Victrex plc. PEEKsil is a trademark of SGE International Pty Ltd Corporation.

All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

Thermo Fisher Scientific Inc. provides this document to its customers with a product purchase to use in the product operation. This document is copyright protected and any reproduction of the whole or any part of this document is strictly prohibited, except with the written authorization of Thermo Fisher Scientific Inc.

The contents of this document are subject to change without notice. All technical information in this document is for reference purposes only. System configurations and specifications in this document supersede all previous information received by the purchaser.

This document is not part of any sales contract between Thermo Fisher Scientific Inc. and a purchaser. This document shall in no way govern or modify any Terms and Conditions of Sale, which Terms and Conditions of Sale shall govern all conflicting information between the two documents.

Release history: Revision A, Apr, 2010; Revision B, Apr 2011; Revision C, Sept 2011; Revision D, Dec 2013

Software version: Thermo Xcalibur 2.1 or later; LC Devices 2.5.0 or later.

**For Research Use Only. Not for use in diagnostic procedures.**

## Regulatory Compliance

Thermo Fisher Scientific performs complete testing and evaluation of its products to ensure full compliance with applicable domestic and international regulations. When the system is delivered to you, it meets all pertinent electromagnetic compatibility (EMC) and safety standards as described in the next section or sections by product name.

Changes that you make to your system may void compliance with one or more of these EMC and safety standards. Changes to your system include replacing a part or adding components, options, or peripherals not specifically authorized and qualified by Thermo Fisher Scientific. To ensure continued compliance with EMC and safety standards, replacement parts and additional components, options, and peripherals must be ordered from Thermo Fisher Scientific or one of its authorized representatives.

### EMC Directive 2004/108/EC

EMC compliance has been evaluated by CTC Analytics for the HTC PAL™ autosampler.

IEC 61326-1:2005

IEC 61326-2-6:2005

EN 61326-1:1997, A1:1998

CISPR 22:2005, A1:2005, A2:2006

FCC Class A, CFR 47 Part 15:2003

### Low Voltage Safety Compliance

Low Voltage Safety Compliance has been evaluated by CTC Analytics for the HTC PAL autosampler.

This device complies with Low Voltage Directive 2006/95/EC and harmonized standard EN 61010-1:2001, IEC 61010-1:2001, ANSI/UL 61010 A-1:2004, CAN/CSA 22.2 61010-1:2004.

## FCC Compliance Statement

THIS DEVICE COMPLIES WITH PART 15 OF THE FCC RULES. OPERATION IS SUBJECT TO THE FOLLOWING TWO CONDITIONS: (1) THIS DEVICE MAY NOT CAUSE HARMFUL INTERFERENCE, AND (2) THIS DEVICE MUST ACCEPT ANY INTERFERENCE RECEIVED, INCLUDING INTERFERENCE THAT MAY CAUSE UNDESIRE OPERATION.



**CAUTION** Read and understand the various precautionary notes, signs, and symbols contained inside this manual pertaining to the safe use and operation of this product before using the device.



## Notice on Lifting and Handling of Thermo Scientific Instruments

For your safety, and in compliance with international regulations, the physical handling of this Thermo Fisher Scientific instrument *requires a team effort* to lift and/or move the instrument. This instrument is too heavy and/or bulky for one person alone to handle safely.

## Notice on the Proper Use of Thermo Scientific Instruments

In compliance with international regulations: This instrument must be used in the manner specified by Thermo Fisher Scientific to ensure protections provided by the instrument are not impaired. Deviations from specified instructions on the proper use of the instrument include changes to the system and part replacement. Accordingly, order replacement parts from Thermo Fisher Scientific or one of its authorized representatives.

## Notice on the Susceptibility to Electromagnetic Transmissions

Your instrument is designed to work in a controlled electromagnetic environment. Do not use radio frequency transmitters, such as mobile phones, in close proximity to the instrument.

For manufacturing location, see the label on the instrument.



## WEEE Compliance

This product complies with the European Union's Waste Electrical & Electronic Equipment (WEEE) Directive 2002/96/EC. It is marked with the following symbol:



Thermo Fisher Scientific has contracted with one or more recycling or disposal companies in each European Union (EU) Member State, and these companies should dispose of or recycle this product. For further information, go to [www.thermoscientific.fr/rohs](http://www.thermoscientific.fr/rohs).

## WEEE Konformität

Dieses Produkt entspricht der EU Waste Electrical & Electronic Equipment (WEEE) Richtlinie 2002/96/EC. Es ist mit dem folgenden Symbol gekennzeichnet:









Thermo Fisher Scientific hat Vereinbarungen mit Verwertungs-/Entsorgungsfirmen in allen EU-Mitgliedsstaaten getroffen, damit dieses Produkt durch diese Firmen wiederverwertet oder entsorgt werden kann. Weitere Informationen finden Sie unter [www.thermoscientific.de/rohs](http://www.thermoscientific.de/rohs).







## Conformité DEEE

Ce produit est conforme avec la directive européenne (2002/96/EC) des Déchets d'Equipements Electriques et Electroniques (DEEE). Il est marqué par le symbole suivant:



Thermo Fisher Scientific s'est associé avec une ou plusieurs sociétés de recyclage dans chaque état membre de l'Union Européenne et ce produit devrait être collecté ou recyclé par celle(s)-ci. Pour davantage d'informations, rendez-vous sur la page [www.thermoscientific.fr/rohs](http://www.thermoscientific.fr/rohs).

CAUTION Symbol	CAUTION	VORSICHT	PRECAUCIÓN	MISE EN GARDE
	<b>Risk electric shock:</b> This instrument uses voltages that can cause electric shock and/or personal injury. Before servicing, shut down the instrument and disconnect it from line power. While operating the instrument, keep covers on. Do not remove the protective covers from the printed circuit board assemblies (PCBAs).	<b>Stromschlaggefahr:</b> Dieses Gerät arbeitet mit Spannungen, die Stromschläge und/oder Personenverletzungen verursachen können. Vor Wartungsarbeiten muss das Gerät abgeschaltet und vom Netz getrennt werden. Betreiben Sie das Gerät nicht mit abgenommenen Abdeckungen. Nehmen Sie die Schutzabdeckungen von Leiterplatten nicht ab.	<b>Riesgo de descargas eléctricas:</b> Este instrumento utiliza voltajes que pueden causar descargas eléctricas y/o lesiones personales. Antes de revisar o reparar el instrumento, apáguelo y desconéctelo de la red eléctrica. Mantenga colocadas las cubiertas mientras se utiliza el instrumento. No retire las cubiertas protectoras del circuito impreso completo (PCBA).	<b>Risque de choc électrique :</b> l'instrument utilise des tensions susceptibles de provoquer une électrocution et/ou des blessures corporelles. Il doit être arrêté et débranché de la source de courant avant toute intervention. Ne pas utiliser l'instrument sans ses couvercles. Ne pas enlever les capots de protection des cartes à circuit imprimé (PCBA).
	<b>Chemical hazard:</b> Wear gloves and other protective equipment, as appropriate, when handling toxic, carcinogenic, mutagenic, corrosive, or irritant chemicals. Use approved containers and proper procedures to dispose of waste oil and when handling wetted parts of the instrument.	<b>Gefahr durch Chemikalien:</b> Tragen Sie beim Umgang mit toxischen, karzinogenen, mutagenen, ätzenden oder reizenden Chemikalien Schutzhandschuhe und weitere geeignete Schutzausrüstung. Verwenden Sie bei der Entsorgung von verbrauchtem Öl und beim Umgang mit medienberührenden Komponenten die vorgeschriebenen Behälter, und wenden Sie ordnungsgemäße Verfahren an.	<b>Peligro por sustancias químicas:</b> Cuando manipule sustancias químicas, tóxicas, carcinogénicas, mutágenas, corrosivas o irritantes, utilice guantes y otro equipo de protección. Utilice siempre recipientes homologados y siga los procedimientos adecuados cuando deseche aceite residual o manipule partes mojadas del instrumento.	<b>Danger lié aux produits chimiques :</b> porter des gants et d'autres équipements de protection appropriés pour manipuler les produits chimiques toxiques, cancérogènes, mutagènes, corrosifs ou irritants. Utiliser des récipients homologués et des procédures adéquates pour la mise au rebut des huiles usagées et lors de la manipulation des pièces de l'instrument en contact avec l'eau.
	<b>Hot surface:</b> Before touching, allow any heated components to cool.	<b>Heiße Oberflächen:</b> Lassen Sie heiße Komponenten vor der Berührung abkühlen.	<b>Superficies calientes:</b> Antes de tocar los componentes calientes, espere a que se enfríen.	<b>Surface chaude :</b> laisser refroidir les composants chauffés avant toute manipulation.
	<b>Flammable substances hazard:</b> Use care when operating the system in the presence of flammable substances.	<b>Gefahr durch entzündbare Substanzen:</b> Beachten Sie die einschlägigen Vorsichtsmaßnahmen, wenn Sie das System in Gegenwart von entzündbaren Substanzen betreiben.	<b>Peligro por sustancias inflamables:</b> Tenga mucho cuidado cuando utilice el sistema cerca de sustancias inflamables.	<b>Danger lié aux substances inflammables :</b> agir avec précaution lors de l'utilisation du système en présence de substances inflammables.
	<b>Risk of eye injury:</b> Eye injury could occur from splattered chemicals, airborne particles, or sharp objects. (Sharp objects that customers might install in the instrument include fused-silica tubing, the autosampler needle, and so on.) Wear safety glasses when handling chemicals or servicing the instrument.	<b>Augenverletzungsrisiko:</b> Verspritzte Chemikalien, Schwebstoffpartikel oder scharfe Objekte können Augenverletzungen verursachen. (Scharfe Objekte, die Kunden möglicherweise im Gerät installieren, sind z. B. Quarzglas-Kapillaren, die Nadel des Autosamplers, usw.) Tragen Sie beim Umgang mit Chemikalien oder bei der Wartung des Gerätes eine Schutzbrille.	<b>Riesgo de lesiones oculares:</b> Las salpicaduras de sustancias químicas, las partículas flotantes en el aire y los objetos afilados pueden causar lesiones oculares. (Entre los objetos afilados que los clientes pueden instalar en el instrumento se encuentran tubos de sílice fundida, agujas del muestreador automático, etc.). Para manipular sustancias químicas o realizar tareas de mantenimiento, utilice gafas de seguridad.	<b>Risque de lésion oculaire :</b> les projections chimiques, les particules en suspension dans l'air et les objets tranchants peuvent entraîner des lésions oculaires. (Les objets tranchants pouvant être installés par les clients dans l'instrument comprennent les tubes en silice fondue, les aiguilles du passeur automatique, etc.). Porter des lunettes de protection lors de toute manipulation de produit chimique ou intervention sur l'instrument.
	<b>General hazard:</b> A hazard is present that is not included in the other categories. This symbol also appears on the instrument. For details about the hazard, refer to the instrument manual. When the safety of a procedure is questionable, contact Technical Support for Thermo Scientific San Jose products.	<b>Allgemeine Gefahr:</b> Es besteht eine weitere Gefahr, die nicht in den vorstehenden Kategorien beschrieben ist. Dieses Symbol wird auch auf dem Gerät angebracht. Einzelheiten zu dieser Gefahr finden Sie in den Gerätehandbüchern. Wenn Sie sich über die Sicherheit eines Verfahrens im Unklaren sind, setzen Sie sich, bevor Sie fortfahren, mit dem technischen Support für Thermo Scientific San Jose Produkte in Verbindung.	<b>Peligro general:</b> Existen peligros que no se incluyen en las otras categorías. Este símbolo también aparece en el instrumento. Si desea obtener más información sobre estos peligros, consulte el manual del instrumento. En caso de duda sobre la seguridad de un procedimiento, póngase en contacto con el personal de servicio técnico de los productos Thermo Scientific San Jose.	<b>Danger d'ordre général :</b> indique la présence d'un risque n'appartenant pas aux catégories citées plus haut. Ce symbole figure également sur l'instrument. Pour plus de détails sur ce danger potentiel, se reporter au manuel de l'instrument. Si la sûreté d'une procédure est incertaine, contacter l'assistance technique pour les produits Thermo Scientific San Jose.

CAUTION Symbol	CAUTION	VORSICHT	PRECAUCIÓN	MISE EN GARDE
	<b>Laser hazard:</b> This instrument uses a laser that is capable of causing personal injury. This symbol also appears on the instrument. For details about the hazard, refer to the instrument manual.	<b>Gefahr durch Laserstrahlen:</b> Der in diesem Gerät verwendete Laser kann zu Verletzungen führen. Dieses Symbol wird auch auf dem Gerät angebracht. Einzelheiten zu dieser Gefahr finden Sie in den Gerätehandbüchern.	<b>Peligro por láser:</b> Este instrumento utiliza un láser que puede producir lesiones personales. Este símbolo también aparece en el instrumento. Si desea obtener más información sobre el peligro, consulte el manual del instrumento.	<b>Danger lié au laser :</b> l'instrument utilise un laser susceptible de provoquer des blessures corporelles. Ce symbole figure également sur l'instrument. Pour plus de détails sur ce danger potentiel, se reporter au manuel de l'instrument.
	<b>Ultra violet light hazard:</b> Do not look directly at the ultra-violet (UV) light or into the UV source. Exposure can cause eye damage. Wear UV eye protection.	<b>Gefahr durch UV-Licht:</b> Richten Sie Ihren Blick nicht direkt auf ultraviolettes Licht (UV-Licht) oder in die UV-Quelle. Dies kann zu Augenschäden führen. Tragen Sie eine UV-Schutzbrille.	<b>Peligro por luz ultravioleta:</b> No mire directamente a una luz ultravioleta (UV) ni a una fuente UV. La exposición puede causar daños oculares. Lleve protección ocular para UV.	<b>Danger lié aux rayons ultraviolets :</b> ne jamais regarder directement la lumière ultraviolette (UV) ou la source d'UV. Une exposition peut entraîner des lésions oculaires. Porter des protections oculaires anti-UV.
	<b>Sharp object:</b> Avoid physical contact with the object.	<b>Scharfes Objekt:</b> Vermeiden Sie den physischen Kontakt mit dem Objekt.	<b>Objeto puntiagudo:</b> Evite el contacto físico con el objeto.	<b>Objet tranchant :</b> éviter tout contact physique avec l'objet.
	<b>Pinch point:</b> Keep hands away from this area.	<b>Quetschgefahr:</b> Halten Sie Ihre Hände von diesem Bereich fern.	<b>Puntos de pincamiento:</b> Mantenga las manos apartadas de esta área.	<b>Risque de pincement :</b> éloigner les mains de cette zone.
	<b>Heavy objects:</b> Never lift or move the instrument by yourself; you can suffer personal injury or damage the equipment. For specific lifting instructions, refer to the instrument manual.	<b>Schweres Objekt:</b> Bewegen und heben Sie das Gerät niemals allein an; dies kann zu Verletzungen oder zur Beschädigung des Geräts führen. Spezifische Anweisungen zum Anheben finden Sie im Gerätehandbuch.	<b>Objeto pesado:</b> Nunca levante ni mueva el instrumento por su cuenta, podría sufrir lesiones personales o dañar el equipo. Para obtener instrucciones específicas sobre levantamiento, consulte el manual del instrumento.	<b>Objet lourd :</b> ne jamais soulever ou déplacer l'instrument seul sous peine de blessure corporelle ou d'endommagement de l'instrument. Pour obtenir des instructions de levage spécifiques, se reporter au manuel de l'instrument.
	<b>Trip obstacle:</b> Be aware of cords, hoses, or other objects located on the floor.	<b>Stolpergefahr:</b> Achten Sie auf Kabel, Schläuche und andere Objekte auf dem Fußboden.	<b>Tropiezo con obstáculos:</b> Tenga en cuenta los cables, mangueras u otros objetos colocados en el suelo.	<b>Risque de trébuchement :</b> faire attention aux câbles, tuyaux et autres objets situés sur le sol.
	When the safety of a procedure is questionable, contact Technical Support for Thermo Scientific San Jose products.	Wenn Sie sich über die Sicherheit eines Verfahrens im unklaren sind, setzen Sie sich, bevor Sie fortfahren, mit Ihrer lokalen technischen Unterstützungsorganisation für Thermo Scientific San Jose Produkte in Verbindung.	En caso de duda sobre la seguridad de un procedimiento, póngase en contacto con el personal de servicio técnico de los productos Thermo Scientific San Jose.	Si la sûreté d'une procédure est incertaine, contacter l'assistance technique pour les produits Thermo Scientific San Jose.

CAUTION Symbol	CAUTION	警告	危険警告
	<p><b>Risk electric shock:</b> This instrument uses voltages that can cause electric shock and/or personal injury. Before servicing, shut down the instrument and disconnect it from line power. While operating the instrument, keep covers on. Do not remove the protective covers from the printed circuit board assemblies (PCBAs).</p>	<p><b>感電の危険性:</b> この機器では、感電および/または身体傷害を引き起こすおそれのある電圧を使用しています。整備点検の前には、機器の電源を切り、電源コードを抜いてください。機器の作動中は、カバーを付けたままにしてください。プリント基板アセンブリ (PCBA) から保護カバーを取り外さないでください。</p>	<p><b>触电危险:</b> 本仪器所用电压可能导致电击或人身伤害。进行维修服务前, 务必关闭仪器电源并断开其电源连接。操作此仪器时, 不要卸下顶盖。勿卸下印刷电路板组件 (PCBA) 的保护盖。</p>
	<p><b>Chemical hazard:</b> Wear gloves and other protective equipment, as appropriate, when handling toxic, carcinogenic, mutagenic, corrosive, or irritant chemicals. Use approved containers and proper procedures to dispose of waste oil and when handling wetted parts of the instrument.</p>	<p><b>化学的危険性:</b> 毒性、発癌性、変異原性、腐食性、または刺激性のある化学薬品を取り扱うときは、必要に応じて手袋などの保護具を着用します。廃油を処分したり、機器の接液部品を取り扱うときは、認可された容器を使用し、適切な手順に従います。</p>	<p><b>化学品危险:</b> 当处理毒性、致癌性、致突变性、腐蚀性或者刺激性化学品时, 佩戴手套和其他保护性设备。当处理浸湿的仪器部件以及废油时, 使用认可的容器和合适的步骤。</p>
	<p><b>Hot surface:</b> Before touching, allow any heated components to cool.</p>	<p><b>高温面:</b> 触れる前に、加熱した部品を冷ましてください。</p>	<p><b>热表面:</b> 待高温部件冷却之后再行维修。</p>
	<p><b>Flammable substances hazard:</b> Use care when operating the system in the presence of flammable substances.</p>	<p><b>可燃性物質の危険性:</b> 可燃性物質があるところでシステムを作動させる場合は十分注意してください。</p>	<p><b>易燃物危险:</b> 在有易燃物质的场地操作该系统时, 务必小心谨慎。</p>
	<p><b>Risk of eye injury:</b> Eye injury could occur from splattered chemicals, airborne particles, or sharp objects. (Sharp objects that customers might install in the instrument include fused-silica tubing, the autosampler needle, and so on.) Wear safety glasses when handling chemicals or servicing the instrument.</p>	<p><b>眼外傷の危険性:</b> 飛散した化学薬品、浮遊粒子、または鋭利な物体によって眼外傷を負うおそれがあります (機器に取り付けられる可能性がある鋭利な物体は、ヒューズドシリカ、オートサンプラーニードルなどです)。化学薬品を取り扱ったり、機器を整備点検するときは、保護メガネを着用します。</p>	<p><b>眼睛伤害风险:</b> 眼睛受伤可能源自飞溅的化学品、空气中的颗粒, 或者锋利的物体。(安装在仪器内的锋利物体包括熔融石英管、自动进样器的进样针等。) 处理化学品或对仪器进行维修服务时, 务必戴上防护眼镜。</p>
	<p><b>General hazard:</b> A hazard is present that is not included in the other categories. This symbol also appears on the instrument. For details about the hazard, refer to the instrument manual. When the safety of a procedure is questionable, contact Technical Support for Thermo Scientific San Jose products.</p>	<p><b>一般的な危険性:</b> それぞれのカテゴリーに当てはまらない危険があります。この標識記号は機器にも表示されています。この危険の詳細については、機器のマニュアルを参照してください。手順の安全性にご不明な点がある場合は、Thermo Scientific San Jose 製品のテクニカルサポートまでお問い合わせください。</p>	<p><b>普通危険:</b> 未归入其他类别的危险。此符号也会在仪器上出现。有关此危险的详细信息, 参阅适当的仪器手册。若对任何步骤的安全事项有疑问, 联系 Thermo Scientific San Jose 产品的技术支持中心。</p>

CAUTION Symbol	CAUTION	警告	危険警告
	<b>Laser hazard:</b> This instrument uses a laser that is capable of causing personal injury. This symbol also appears on the instrument. For details about the hazard, refer to the instrument manual.	<b>レーザー光線の危険性:</b> この機器では、身体傷害を引き起こすおそれのあるレーザーを使用しています。この標識記号は機器にも表示されています。この危険の詳細については、機器のマニュアルを参照してください。	<b>激光危险:</b> 本仪器所用激光会导致人身伤害。此符号也会在仪器上出现。有关此危险的详细信息，参阅适当的仪器手册。
	<b>Ultra violet light hazard:</b> Do not look directly at the ultra-violet (UV) light or into the UV source. Exposure can cause eye damage. Wear UV eye protection.	<b>紫外光の危険性:</b> 紫外 (UV) 光または UV 光源を直接見ないでください。照射によって眼損傷を引き起こすおそれがあります。UV 保護メガネを着用します。	<b>紫外光危険:</b> 不要直视紫外 (UV) 光或者紫外光源。直视可能导致眼睛伤害。佩戴紫外线防护眼镜。
	<b>Sharp object:</b> Avoid physical contact with the object.	<b>鋭利な物体:</b> 物体との身体的接触を避けてください。	<b>锋利物体:</b> 避免直接接触锋利的物体。
	<b>Pinch point:</b> Keep hands away from this area.	<b>ピンチポイント:</b> この部分には手を挟まれないようにしてください。	<b>夹点:</b> 勿将手放在此部位。
	<b>Heavy objects:</b> Never lift or move the instrument by yourself; you can suffer personal injury or damage the equipment. For specific lifting instructions, refer to the instrument manual.	<b>重量物:</b> 1人で機器を持ち上げたり移動しないでください。身体傷害を負ったり、機器を損傷するおそれがあります。具体的な持ち上げ方法については、機器のマニュアルを参照してください。	<b>重物:</b> 切勿独自提起或移动本仪器; 可能遭受人身伤害或损坏仪器。有关具体的提起说明, 参阅仪器手册。
	<b>Trip obstacle:</b> Be aware of cords, hoses, or other objects located on the floor.	<b>作業の障害物:</b> 床にあるコード、ホース、その他の物体に注意してください。	<b>绊倒危険:</b> 注意地面上的线、管或其他物品。
	When the safety of a procedure is questionable, contact Technical Support for Thermo Scientific San Jose products.	手順の安全性にご不明な点がある場合は、Thermo Scientific San Jose 製品のテクニカルサポートまでお問い合わせください。	如对安全程序有疑问, 联系 Thermo Scientific San Jose 产品的技术支持中心。

# Contents

	<b>Preface</b> .....	<b>xv</b>
	Related Documentation .....	xv
	Cautions and Special Notices .....	xvi
	Contacting Us .....	xvii
<b>Chapter 1</b>	<b>Introduction</b> .....	<b>1</b>
	Open Autosampler Components .....	1
	Main Components .....	2
	Dynamic Load and Wash (DLW) .....	3
	Specifications .....	6
	Instrument Dimensions .....	8
	Operating and Environmental Requirements .....	9
	Sound Pressure Level .....	9
<b>Chapter 2</b>	<b>Installing the Accela Open Autosampler</b> .....	<b>11</b>
	Installing the Autosampler and DLW .....	11
	Unpacking the Components .....	12
	Assembling the Accela Open Autosampler .....	14
	Installing the DLW Option .....	25
	Electrical Connections .....	36
	System Synchronization Connections .....	39
	Accela Open Autosampler Object Positions .....	40
	Defining Object Positions .....	40
	Description of Object Positions .....	41
	Injection Valve .....	50
	Valve Drives and Valves – General Remarks .....	50
	Injection Valve Flow Path .....	50
<b>Chapter 3</b>	<b>Operating the Accela Open Autosampler</b> .....	<b>55</b>
	Control Terminal .....	55
	Menu Screens .....	56
	Function Keys .....	57
	ESC and STOP Keys .....	57
	Scroll Knob and Enter Button .....	57



Methods	57
Creating Methods	57
Editing and Viewing Methods	58
Deleting Methods	59
Job and Job Queue	59
Building and Starting a Job Queue	60
Canceling a Job Queue	61
Restarting a Canceled Job Queue	61
Utility Functions	61
Syringe	62
Tray	63
Injector	64
Wash Station	65
Vial	66
Dilutors	66
Tools	67
Logfile	68
Info Functions	68
Hardware	69
Software	69
Maintenance	69
Free Objects/Free Items	70
Setup Functions	71
Sounds	71
Time	71
Objects	72
Communicating with the Computer	72
<b>Chapter 4</b>	<b>Operating the Temperature Controllable Stack Coolers</b>
	<b>75</b>
Temperature Setting, Control, and Alarm	75
Temperature Setting	75
Temperature Control	76
Temperature Alarm	77
Temperature Stability	77
Condensation Build-Up	78
Condensation Build-Up in Accela Open Autosampler Stack Coolers	78
Condensation Build-Up in PAL Tray Coolers	79
<b>Chapter 5</b>	<b>Maintenance</b>
	<b>81</b>
<b>Chapter 6</b>	<b>Troubleshooting</b>
	<b>85</b>
Troubleshooting Using Observed Symptoms or Error Messages	85
Cooled Stack Power Supply Error Messages	90

Troubleshooting Considerations . . . . .	90
Tubing Connections . . . . .	90
Solvent Delivery Pumps Not Priming . . . . .	95
Solvent Delivery Pump Not Functioning . . . . .	95
Standard Chromatographic Tests . . . . .	96
Carryover Test with UV-Detection . . . . .	97
Repeatability Test with UV-Detection. . . . .	98
<b>Chapter 7 Replaceable Parts . . . . .</b>	<b>101</b>
DLW Pump Module . . . . .	102
DLW Wash Station . . . . .	102
DLW Syringe . . . . .	102
DLW Syringe Holder Assembly . . . . .	103
DLW Option Spare Parts Ordering Information. . . . .	103
DLW-2 Option Spare Parts Ordering Information . . . . .	108
Control-xt Board . . . . .	109
Injection Valve and Valve Rotor . . . . .	110
Cleaning or Replacing Injection Valves . . . . .	110
Cleaning or Replacing Valve Rotors . . . . .	112
Injection Unit . . . . .	112
<b>Appendix A Reference Information . . . . .</b>	<b>113</b>
Glossary . . . . .	114
Naming Convention . . . . .	116
Special Functions . . . . .	117
Accessing Extended User Mode . . . . .	117
Section F3-Setup . . . . .	117
Section F3-Setup/Objects . . . . .	124
PAL Firmware Overview . . . . .	136
PAL Firmware Error Codes . . . . .	140
External Connectors for Accela Open Autosampler . . . . .	145
System Limitations . . . . .	152
Accela Open Autosampler . . . . .	152
Sample Volume . . . . .	152
Wetted Parts and Material. . . . .	152
Tubing Internal Diameter (ID) and System Backpressure . . . . .	152
Interfacing the Open Autosampler to Other Devices. . . . .	152
<b>Appendix B PAL Loader . . . . .</b>	<b>155</b>
Introduction . . . . .	155
New Features of PAL Loader Software 2.1.x . . . . .	156
Installation of PAL Loader Software . . . . .	156
Requirements . . . . .	156
PAL Loader Software Version 2.1.x Installation. . . . .	157

Multiple Shortcuts Assignment . . . . .	158
Dedicated Shortcut for Serial Communication . . . . .	159
Creating Multiple Shortcuts . . . . .	160
PAL Loader Software Operation . . . . .	161
Starting the PAL Loader Software . . . . .	161
Setting up COM Ports . . . . .	162
Serial Communication. . . . .	164
Reset PAL-xt to Serial Communication. . . . .	165
Backing Up the PAL Firmware . . . . .	166
Updating the PAL Firmware . . . . .	168
Troubleshooting . . . . .	171
Serial Communication with PAL Loader Software. . . . .	171
Reset PAL to Serial Communication . . . . .	172
Erratic Behavior of Software or Flash Memory . . . . .	176
Failure to Complete BACKUP or UPDATE Process. . . . .	176
Limitations . . . . .	177
<b>Appendix C PAL Object Manager. . . . .</b>	<b>179</b>
Installing PAL Object Manager . . . . .	180
Verifying System Requirements. . . . .	180
Installing the Object Manager Software to the Local Hard Drive . . . . .	180
Installing the PAL Object Lists . . . . .	181
Getting Started with PAL Object Manager . . . . .	181
Setting Up the Communication Mode . . . . .	182
Troubleshooting the Serial Communication Mode . . . . .	182
Using PAL Object Manager to Change Your System Configuration . . . . .	183
<b>Appendix D Column Switching Option . . . . .</b>	<b>187</b>
Installing the Column Switching Valves . . . . .	187
Installing the 6-Port Valve. . . . .	188
Installing the 10-Port Valve. . . . .	189
Installing the Software Tools . . . . .	191
Backing Up the System . . . . .	192
Installing the Column Switch Option . . . . .	193
Installing a New Backup File to the Autosampler . . . . .	193
Installing Special Objects to the Autosampler . . . . .	194
Verifying the Valve Configuration. . . . .	196
Installing the Macros and Methods . . . . .	197
Creating a Method in Xcalibur . . . . .	198
<b>Appendix E Changing the DLW Holding Loop . . . . .</b>	<b>201</b>
<b>Index . . . . .</b>	<b>209</b>

# Preface

This manual describes the features, installation, and maintenance of the Accela™ Open Autosampler.

## Contents

- [Related Documentation](#)
- [Cautions and Special Notices](#)
- [Contacting Us](#)

### ❖ To suggest changes to the documentation or to the Help

Complete a brief survey about this document by clicking the button below. Thank you in advance for your help.



## Related Documentation

The Accela Open Autosampler includes complete documentation. In addition to this guide, you can also access the following documents as PDF files from the data system computer:

- *Accela LC System Preinstallation Requirements Guide*
- *Accela UHPLC System Getting Connected Guide*
- *Accela Open Autosampler User Guide for LC Devices*

❖ **To view the product manuals**

From the Microsoft™ Windows™ taskbar, do the following:

- For a Thermo Scientific mass spectrometer, choose **Start > All Programs > Thermo Instruments** and so on.
- For an LC instrument controlled by a Thermo software application, choose **Start > All Programs > Thermo Instruments > Manuals > LC Devices > Accela** and so on.

For access to the application Help, follow this procedure.

❖ **To view application-specific Help**

- From the Accela Open Autosampler window, choose **Help > Accela Open Autosampler Help**.
- If information about setting parameters is available for a specific view, page, or dialog box, click **Help** or press the F1 key for information about setting parameters.

For more information, visit [www.thermoscientific.com](http://www.thermoscientific.com).

## Cautions and Special Notices

Make sure you follow the cautions and special notices presented in this guide. Cautions and special notices appear in boxes; those concerning safety or possible system damage also have corresponding caution symbols.

This guide uses the following types of cautions and special notices.



**CAUTION** Highlights hazards to humans, property, or the environment. Each CAUTION notice is accompanied by an appropriate CAUTION symbol.

**IMPORTANT** Highlights information necessary to prevent damage to software, loss of data, or invalid test results; or might contain information that is critical for optimal performance of the system.

**Note** Highlights information of general interest.

**Tip** Highlights helpful information that can make a task easier.

## Contacting Us

There are several ways to contact Thermo Fisher Scientific for the information you need.

### ❖ To contact Technical Support

Phone	800-532-4752
Fax	561-688-8736
E-mail	<a href="mailto:us.techsupport.analyze@thermofisher.com">us.techsupport.analyze@thermofisher.com</a>

### ❖ To contact Customer Service for ordering information

Phone	800-532-4752
Fax	561-688-8731
E-mail	<a href="mailto:us.customer-support.analyze@thermofisher.com">us.customer-support.analyze@thermofisher.com</a>
Web site	<a href="http://www.thermoscientific.com/en/support-landing/support.html">http://www.thermoscientific.com/en/support-landing/support.html</a>

### ❖ To copy manuals from the Internet

Go to [mssupport.thermo.com](http://mssupport.thermo.com), agree to the Terms and Conditions, and then click **Customer Manuals** in the left margin of the window.

### ❖ To suggest changes to the Help

- Fill out a reader survey online at [www.surveymonkey.com/s/PQM6P62](http://www.surveymonkey.com/s/PQM6P62).
- Send an e-mail message to the Technical Publications Editor at [techpubs-lcms@thermofisher.com](mailto:techpubs-lcms@thermofisher.com).





# Introduction

For an overview of the Accela Open Autosampler, read the following topics.

## Contents

- [Open Autosampler Components](#)
- [Specifications](#)
- [Instrument Dimensions](#)
- [Operating and Environmental Requirements](#)
- [Sound Pressure Level](#)

## Open Autosampler Components

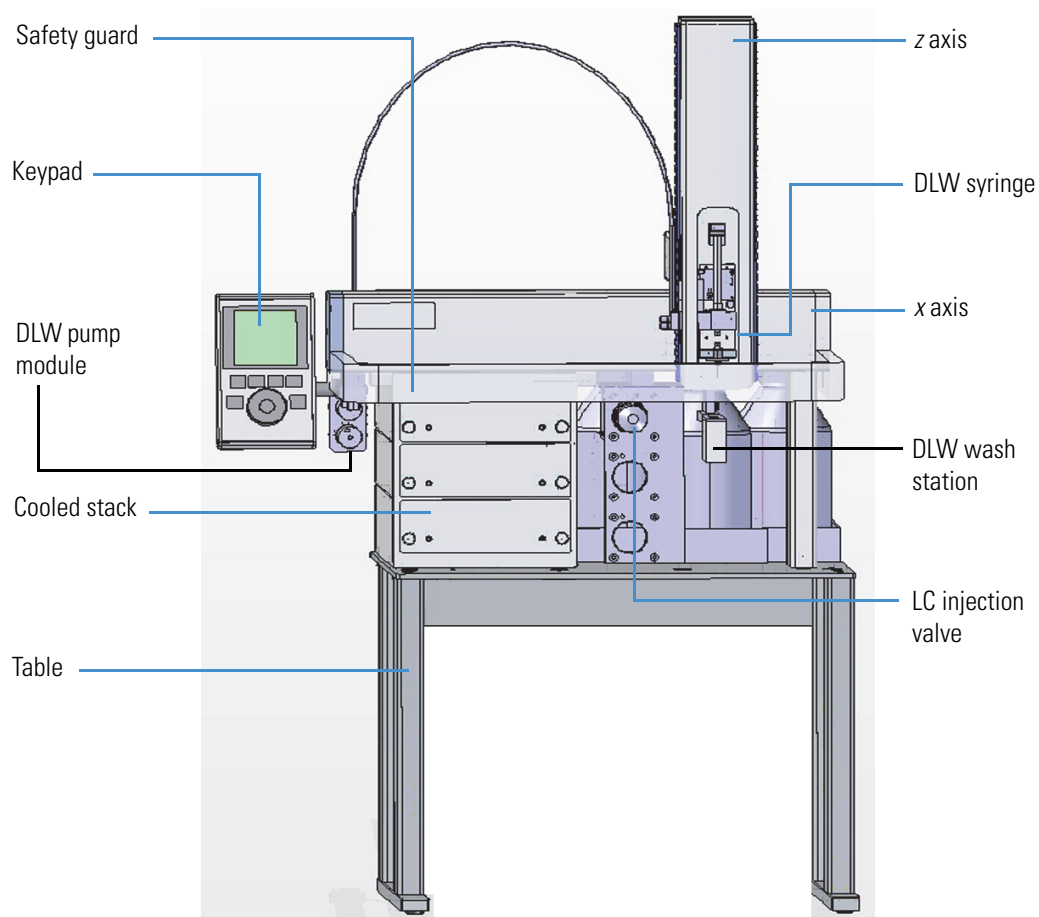
This section introduces you to the major components of the Accela Open Autosampler.

- [Main Components](#)
- [Dynamic Load and Wash \(DLW\)](#)

## Main Components

Figure 1 shows the location for the major system components for the Accela Open Autosampler.

**Figure 1.** Accela Open Autosampler major system components



The main components of the Accela Open Autosampler are as follows:

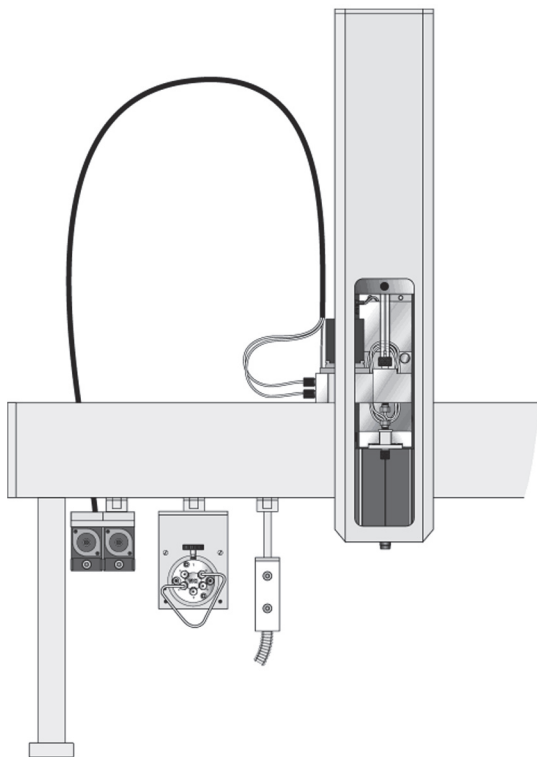
- $x$ -,  $y$ -axes assembly
- Injection unit,  $z$  axis
- Syringe kit for liquid injections
- Fast wash station for two different solvents
- Keypad terminal
- Safety guard
- Standard stand-alone supports, 215 mm

To set up the Accela Open Autosampler system, you need a tray set (tray holder and tray) and a stack. You can choose from many optional types of tray sets. For instructions, refer to the *Accela Open Autosampler User Guide*.

## Dynamic Load and Wash (DLW)

This section contains an overview of the Dynamic Load and Wash (DLW) option (Figure 2).

**Figure 2.** Dynamic Load and Wash (DLW)



The DLW option represents a new wash station concept that combines an injection cycle with wash steps. The linked combination of the two steps, which are usually separate, minimizes cycle time and carryover.

There are two characterizing features for the DLW option:

- The sample solution never contacts the syringe itself; it is held between the holding loop.
- Wash solvents are pumped from back to front into the DLW system to intensely flush all critical parts that are in contact with the sample.

The DLW option consists of two self-priming micro pumps (mounted on a dedicated bracket), and the wetted parts are Ryton™ PPS and Kalrez™ (FFPM). The pump IN ports are connected to the wash solvent bottles and the OUT ports are connected to the DLW manifold, which is part of the assembly of the dedicated DLW syringe holder. A holding loop separates the syringe and the DLW actuator to avoid sample from contacting these parts. For a detailed list of the DLW components, see “DLW-2 components” on page 14.

The syringe and holding loop are preloaded with wash solvent #1 at the start. The sample is picked up and remains separated from wash solvent #1 by an air gap. After loading the loop and injection, wash solvent #1 is pushed into the system, followed directly by wash solvent #2 to flush the critical valve paths.

The DLW syringe assembly is moved to the wash station for further cleaning steps and for preparing the Syringe and holding loop for the next cycle.

There has been continuous development on the DLW option. The current version, DLW-2, has the following improvements over the DLW option:

- Holding Loop

The DLW-2 Holding Loop consists of one single piece from the needle to the loop, which eliminates several tube connections and replaces the needle and flow diverter. The Holding Loop is made of high-quality stainless steel and the inner surface is passivated with acid.

- DLW Syringe Holder back plate

The DLW-2 Syringe Holder back plate has changed in size and form to allow a replacement of the Holding Loop. The Actuator is supported from the back side to provide higher mechanical stability.

- DLW Pump Assembly

The DLW-2 Pump Assembly is equipped with a housing to protect the electrical parts from solvent splashes.

**Note** For a list of the DLW-2 spare parts, see [“DLW-2 Option Spare Parts Ordering Information”](#) on page 108.

## Differences Between DLW and DLW-2

Table 1 shows the differences between the critical parts or parameters of the DLW and DLW-2 options.

**Table 1.** Differences Between DLW and DLW-2

Part or parameter	DLW option	DLW-2 option
<b>Holding Loop</b>		
ID	0.50 mm	0.41 mm
Length	550 mm	855 mm
Material	FEP tubing	Stainless steel, passivated
OD	1/16 in. (1.58 mm)	0.72 mm (Gauge 22)
Volume	108 µL	118 µL
<b>Syringe Needle</b>		
Length	51 mm	Not applicable, integrated in DLW-2 Holding Loop
Gauge	22	
Material	Stainless steel	
<b>DLW Flow Diverter</b>		
Length	20.5 mm	Not applicable, integrated in DLW-2 Holding Loop
OD	0.72 mm (Gauge 22)	
ID	0.41 mm	
Material	Stainless steel	
<b>DLW Internal Volume</b>		
Total delay volume	205 µL + Injection loop volume (204.7)	208 µL + Injection loop volume
	Manifold: 90 µL	Manifold: 90 µL
	Holding Loop: 108 µL	Holding Loop: 118 µL
	Syringe Needle: 6.7 µL	
	Installed Injection Loop	Installed Injection Loop

## Specifications

Table 2 lists the specifications for the Accela Open Autosampler.

**Table 2.** Accela Open Autosampler specifications (Sheet 1 of 3)

Component	Specification
Sample capacity	<ul style="list-style-type: none"> <li>• Six trays with 54 positions for 2 mL vials</li> <li>• Six deepwell or standard microtiter-plates with 96 positions</li> <li>• Six microtiter-plates with 384 positions</li> <li>• Six trays with 15 positions for 10 mL vials</li> </ul>
Thermostatted sample tray	<ul style="list-style-type: none"> <li>• 4 to +40 °C in 0.1 °C increments Peltier element</li> <li>• Temperature cooling: The specification for the stack cooler DW temperature cooling is defined as: <ul style="list-style-type: none"> <li>– <math>T\Delta = \geq 14.0</math> °C within 60 minutes</li> <li>– <math>T\Delta = T_{\text{Roomtemperature}} - T_{\text{Stack Cooler DW}}</math></li> </ul> </li> </ul> <p>You must reach a temperature of 14.0 °C between room temperature (22 ±2 °C) and the stack cooler DW within 60 minutes.</p> <p>The control unit and the resulting temperature display on the stack cooler DW (display) are determinative; no comparison to an independent temperature measurement device is necessary.</p>
Syringe sizes	<ul style="list-style-type: none"> <li>• 10, 25, 50, 100, 250, 500, 1000, 2500, and 5000 µL</li> <li>• Standard DLW syringe: 100 µL</li> </ul>
Liquid injection range	<ul style="list-style-type: none"> <li>• 10–100 µL with standard 100 µL syringe</li> <li>• Down to 0.1 µL with solvent sandwich technique</li> <li>• Down to 0.5 µL with optional 4-port internal loop valve</li> <li>• Up to 5000 µL with optional larger syringe and loop</li> </ul>
Replicate injections	<ul style="list-style-type: none"> <li>• 1–99 from one vial</li> </ul>
Minimum sample volume	<ul style="list-style-type: none"> <li>• 1 µL from a 3 µL sample in 1 mL tapered micro vial</li> </ul>
LC injector	<ul style="list-style-type: none"> <li>• Electrically actuated fast switching valve drive</li> <li>• 6-port Injection valve</li> <li>• VICI™/Valco™ Cheminert™ (P/N 00950-01-00336)</li> </ul> <p>Standard valve and other valves are available upon request.</p>

**Table 2.** Accela Open Autosampler specifications (Sheet 2 of 3)

Component	Specification
Wash station	<p>Fast wash station with two different wash DLW Pumps:</p> <ul style="list-style-type: none"> <li>• Two pieces. solenoid diaphragm pump</li> <li>• Self-priming               <ul style="list-style-type: none"> <li>– Flow rate range: 100 mL/min at atmospheric pressure (Pump itself, no connections to HPLC system)</li> <li>– Wetted parts: Pumps: Ryton PPS, and Kalrez (FFPM)</li> </ul> </li> <li>• Transfer tubing kit: Two pieces. PFA tubes 1/8 in. including connection fittings</li> <li>• Solvent Reservoir: 1000 mL borosilicate glass including 10 µm solvent filter</li> </ul>
Wash station	<ul style="list-style-type: none"> <li>• DLW Wash Station:               <ul style="list-style-type: none"> <li>– Wash station with two solvent ports and one waste port.</li> <li>– Two additional horizontal waste inlet ports in front to be combined with valve waste port.</li> </ul> </li> <li>• Wetted parts:               <ul style="list-style-type: none"> <li>– Wash station block: PVDF; inserts: sst 1.4404</li> <li>– Connections at bottom of inserts: PEEK™</li> </ul> </li> </ul> <p>Waste tube: Polyethylene (PE)</p>



**Table 2.** Accela Open Autosampler specifications (Sheet 3 of 3)

Component	Specification
Wash station	<ul style="list-style-type: none"> <li>• DLW Syringe assembly:               <ul style="list-style-type: none"> <li>– Syringe: Gastight Syringe 100 µL with Valflon tip Wetted parts: Glass, PTFE and Valflon</li> <li>– Syringe needle: Gauge 22 (18/25 mm OD/0.41 mm ID) Length: 51 mm, PTFE seal</li> </ul> </li> <li>• DLW Actuator/solenoid: 2/2-way flipper solenoid valve               <ul style="list-style-type: none"> <li>– Wetted parts: Body: PEEK Seal material: FFKM (Simriz™)</li> </ul> </li> </ul>
Wash station	<ul style="list-style-type: none"> <li>• DLW Manifold Solvent Selector module and Syringe Adapter assembly:               <ul style="list-style-type: none"> <li>– Wetted parts: Block: sst 1.4435 Perfluor (O-ring)</li> </ul> </li> <li>• Holding Loop: FEP Tubing OD 1/16 in., ID 0.50 mm, length 55 cm Loop Volume: 108 µL</li> <li>• DLW Flow Diverter               <ul style="list-style-type: none"> <li>– Tube dimensions: Gauge: 22 (OD 18/25 mm/ID 0.41 mm) Length: 23 mm Wetted part: Stainless chromium steel</li> </ul> </li> <li>• DLW Needle Holder assembly Wetted part: sst 1.4435</li> <li>• Standard Configuration Accela 1 (1250 bar), Accela 2 (1250 bar with 6-drawer cooled stack), Accela 3 (QuickQuan™), Accela 4 (EQuan™)</li> </ul>

## Instrument Dimensions

The dimensions of the Accela Open Autosampler are as follows:

- Height: 1045 mm (41.1 in.)
- Width: 676 mm (26.6 in.)
- Depth: 717.5 mm (28.3 in.)
- Weight: 38.5 kg (84.9 lb)

## Operating and Environmental Requirements

Table 3 lists the operating and environmental requirements of the Accela Open Autosampler.

**Table 3.** Operating and environmental requirements

Parameter	Requirements
Operating temperature range	4 –40 °C (39–104 °F)
Maximum relative humidity	75%, non-condensing
Bench space	<ul style="list-style-type: none"> <li>• At least 24 cm (10 in.) at the back</li> <li>• Access to power switches and power cords</li> <li>• Clean, level and smooth surface</li> <li>• Solid bench plate</li> </ul>
Vibration	Negligible
Static electricity	Negligible

## Sound Pressure Level

Table 4 lists the sound pressure level of the Accela Open Autosampler.

**Table 4.** Sound pressure level

Parameter	Requirements
Sound pressure level	<p>Measured value: 62 dBA (PAL System used for measurement)</p> <p>One meter from the equipment in the direction of maximum sound pressure level</p> <p>According to standard UL 610107A-1, 1st edition, clause 12.5 Limit &lt; 85 dBA</p> <p>dBA = A weighted sound pressure level</p>

**Note** Thermo Fisher Scientific reserves the right to make improvements, changes, or both to the product specifications without notice.



# Installing the Accela Open Autosampler

This chapter describes the installation, contact closure cable connection, computer connection, object positions, and injection valve of the Accela Open Autosampler.

## Contents

- [Installing the Autosampler and DLW](#)
- [System Synchronization Connections](#)
- [Accela Open Autosampler Object Positions](#)
- [Injection Valve](#)

## Installing the Autosampler and DLW

These topics describe how to unpack the components and assemble the open autosampler, install the dynamic load and wash hardware (DLW), electrical connections, and the cool stack.

- [Unpacking the Components](#)
- [Assembling the Accela Open Autosampler](#)
- [Installing the DLW Option](#)
- [Electrical Connections](#)

### Unpacking the Components

An Accela Open Autosampler system is shipped in three boxes. The contents of these boxes are as follows:

Box 1: Contains the  $x$ -,  $y$ -axes assembly, the injection unit, keypad terminal, stand-alone supports, connecting cables, power supply, syringe kit, wash station assembly, safety guard, and miscellaneous parts. See “Open Autosampler Components” on [page 1](#).

Box 2: Contains the DLW injection kit.

Box 3: Contains the cooled stack and the Accela Open Autosampler table.

#### ❖ To unpack the components

1. Open Box 1 and first remove the accessory boxes and the injection unit before attempting to remove the  $x$ -,  $y$ -axes assembly.
2. Carefully lift the  $x$ -,  $y$ -axes assembly and remove it from the box. Hold the  $y$  axis in place while you are removing the assembly from the box. Set the  $x$ -,  $y$ -axes assembly on a bench.
3. Unpack the remaining small boxes and any other accessories.
4. When placing the Open Autosampler onto a stable surface, make sure that no objects interfere with either the  $y$  axis or the injection unit throughout the entire range of potential movement.

The Accela Open Autosampler Dynamic Load and Wash (DLW) option ships in one box (Box 2). Check for the following items:

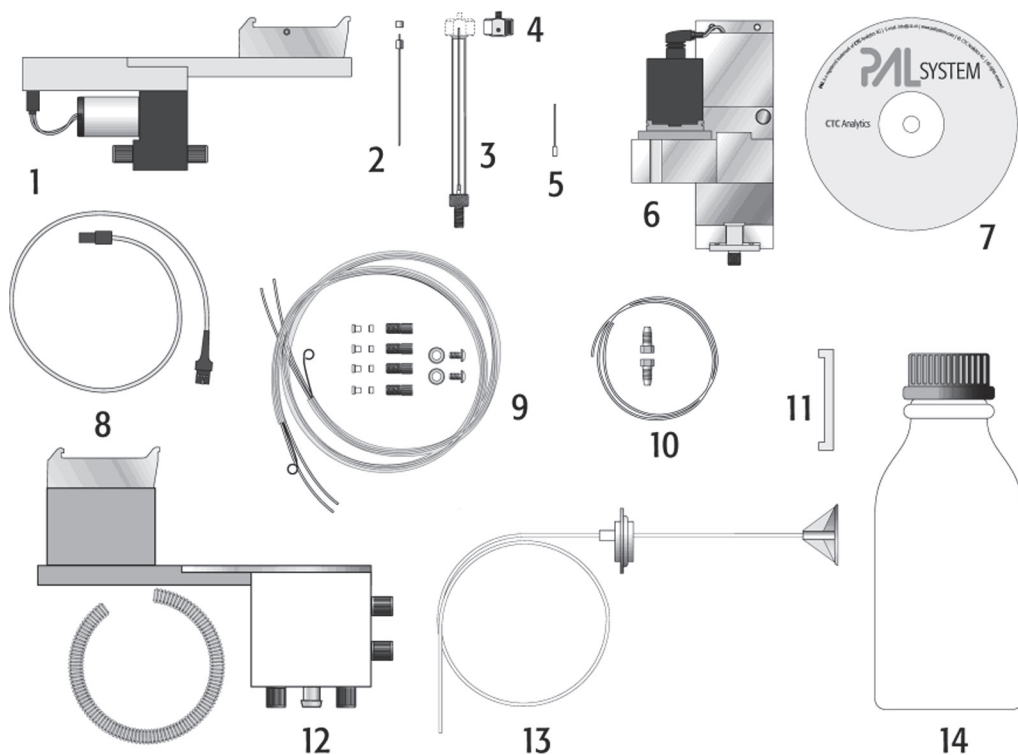
1. DLW Pump Module
2. DLW Syringe Needle
3. DLW Syringe
4. DLW Plunger Holder
5. DLW Flow Diverter (installed in DLW Syringe assembly)
6. DLW Syringe Holder assembly
7. CD-ROM
8. Wash station cable to PAL System
9. DLW Tubing Kit
10. DLW Holding Loop (installed in DLW Syringe assembly)
11. Needle guide length tool
12. DLW Wash Station incl. waste tube

13. 2 pcs. Solvent bottle transfer line including PEEK solvent filter, 10  $\mu\text{m}$
14. 2 pcs. 1 liter solvent reservoir bottles

Figure 3 illustrates the various DLW components as shown in the list above.

For details on how to install the DLW, see “Installing the DLW Option” on page 25. For a list of DLW spare parts and ordering information, see “DLW Option Spare Parts Ordering Information” on page 103.

**Figure 3.** DLW components



The DLW-2 option ships in one box. Check for the following items (Figure 4):

1. DLW Pump Module
2. DLW Syringe
3. DLW Plunger Holder
4. DLW Syringe Holder assembly
5. CD-ROM
6. Wash station cable to PAL System
7. DLW Tubing Kit
8. DLW Holding Loop (installed in DLW Syringe assembly)

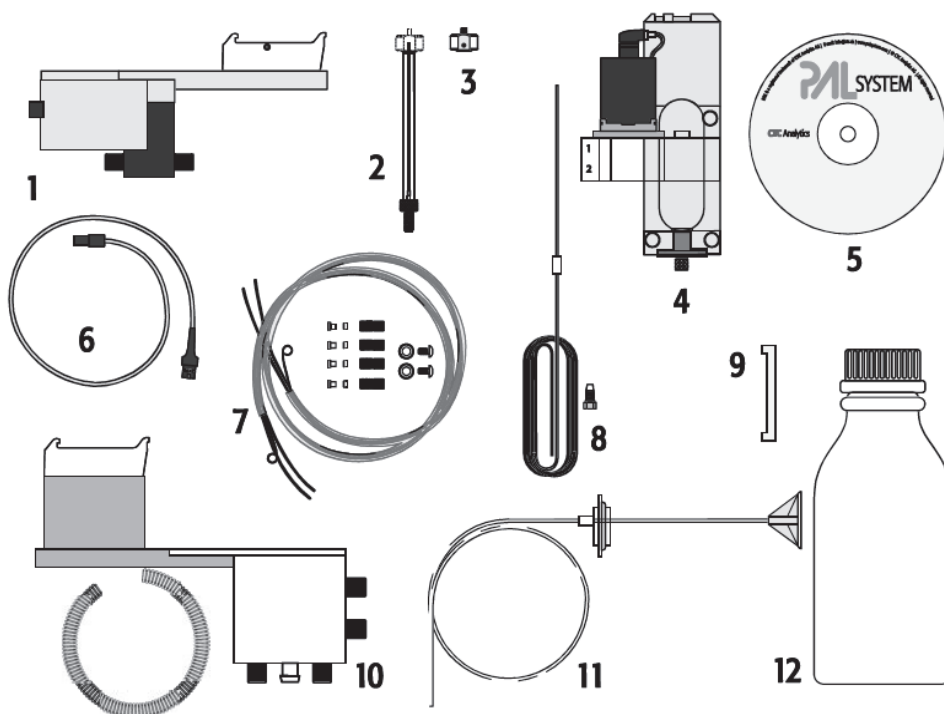
## 2 Installing the Accela Open Autosampler

Installing the Autosampler and DLW

9. Needle guide length tool
10. DLW Wash Station incl. waste tube
11. 2 pcs. Solvent bottle transfer line including glass filter, solvent inlet, 40  $\mu\text{m}$  pore size
12. 2 pcs. 1 liter solvent reservoir bottles

For a list of DLW-2 spare parts and ordering information, see “DLW-2 Option Spare Parts Ordering Information” on page 108.

**Figure 4.** DLW-2 components



## Assembling the Accela Open Autosampler

To assemble the Accela Open Autosampler, follow these procedures:

- [Installing the Cool Stack](#)
- [Installing the Injection Unit](#)
- [Installing the Keypad Terminal](#)
- [Installing the Power Supply](#)



- Installing the LC Injection Valve
- Mounting the DLW Pump Holder

**Tip** Use the screwdrivers in Box 1 to install the different parts with the supplied screws.

## Installing the Cool Stack

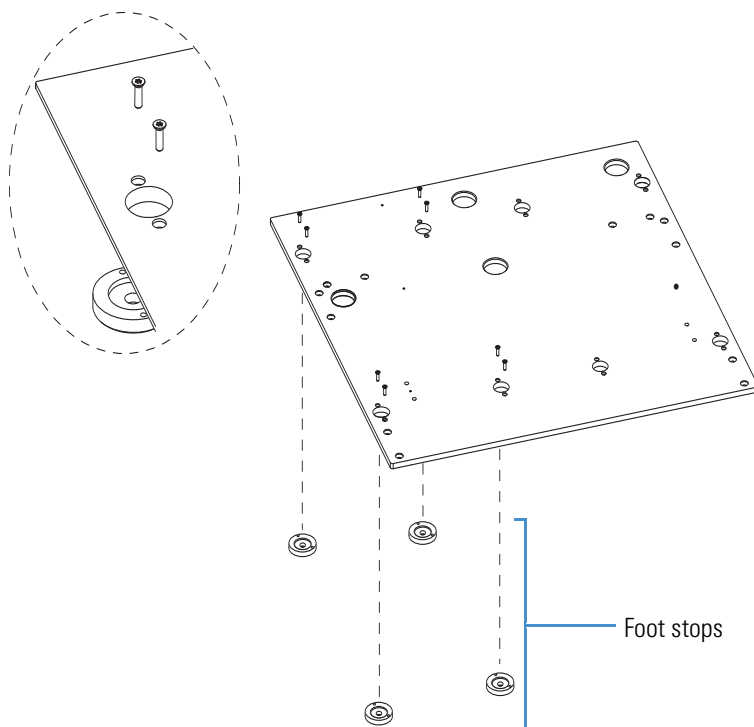
Follow these instructions to install the cool stack on the table and the  $x$  axis on the cool stack.

**Note** A standard table plate is shipped with a left-side cool stack installation. This configuration works well with the TSQ Quantum™, Vantage™, and Velos™ mass spectrometers. On the Exactive™ mass spectrometer, the ion source is mounted to the right side of the mass spectrometer. To minimize the distance between the autosampler injection valve and the Exactive ion source, install the stack cooler on the right side of the table.

### ❖ To install the cool stack on the table

1. Decide on which side of the table you want the cool stack to be located.
2. If you want to relocate the cool stack, remove and relocate the foot stops to the desired location (Figure 5).

**Figure 5.** Relocating the foot stops



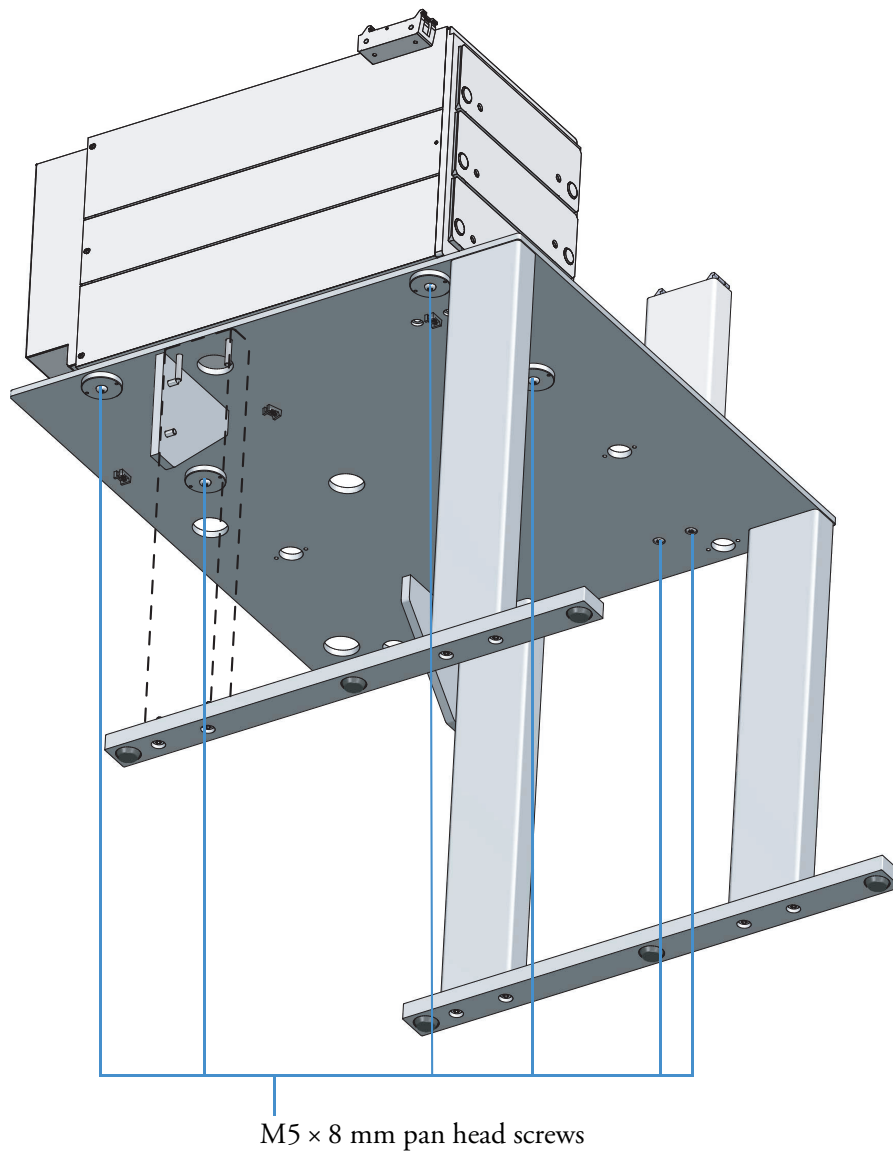
3. Place the cool stack onto the table in the desired location, ensuring that the mounting feet reside within the designated cutouts resting on the foot stops in the table.

## 2 Installing the Accela Open Autosampler

Installing the Autosampler and DLW

- Using the supplied M5 × 8 mm pan head screws (qty 4), secure the cool stack in place (Figure 6).

**Figure 6.** Installing the cool stack to the table

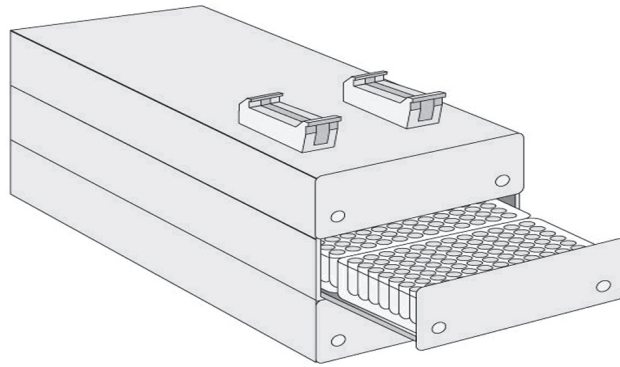


- Using the supplied M5 × 8 mm pan head screws (qty 2), attach the remaining leg for the Open Autosampler itself to the opposite side of the cool stack (Figure 6).

❖ **To install the x axis on the cool stack**

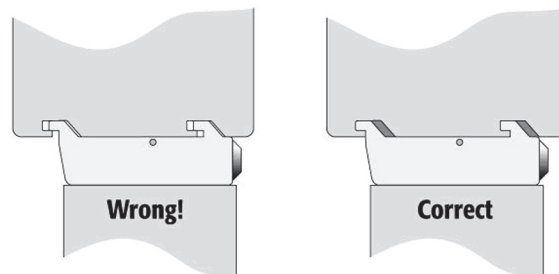
1. Loosen the two Torx™ screws on the three mounting clamps (2 located on top of the stack and 1 on top of the autosampler leg). [Figure 7](#) shows the stack model with 3 drawers.

**Figure 7.** A microplate stack with three drawers



2. Carefully lift the x-axis assembly on top of the stack with the mounting clamp teeth fitting into the grooves on the bottom of the x axis.
3. Ensure the clamps fit completely into the grooves. Alternately tighten the three Torx screws until the three mounting clamps are firmly in place.
4. Ensure the stack clamps are correctly attached to the x axis ([Figure 8](#)). Also ensure the aligned attachment of the x axis to both the cool stack and autosampler leg.

**Figure 8.** Attachment of stack clamps

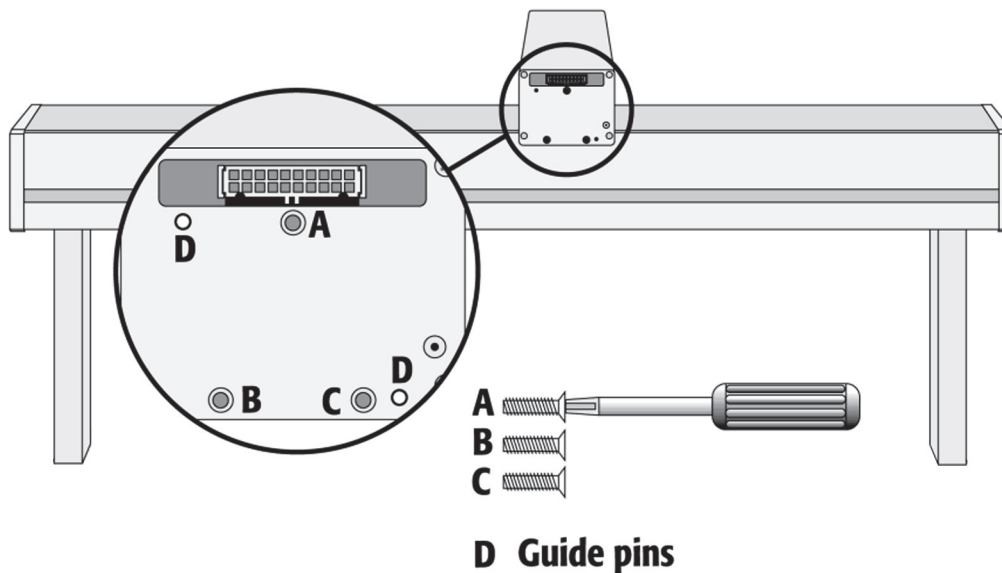


## Installing the Injection Unit

### Tip

- Install the *x*-axis stand before you install the injection unit.
- Install the injection unit with caution. When installing it for the first time, have someone hold it in place while you insert the mounting screws.

**Figure 9.** Attaching the Open Autosampler injection unit

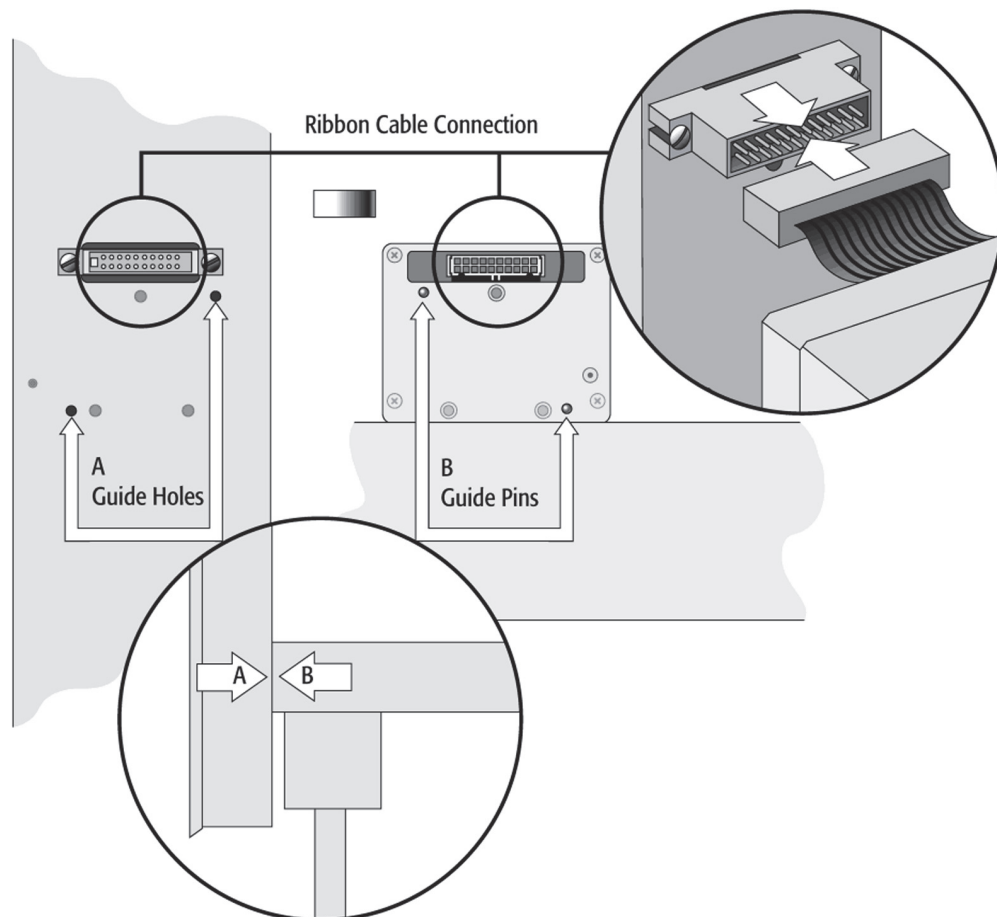


### ❖ To install the injection unit

1. Remove the three Torx mounting screws A, B, and C, used to fix the injection unit to the *y* axis (Figure 9).
2. Connect the ribbon cable A protruding from the front end of the *y* axis to the corresponding connector on the injection unit (Figure 10).

**Tip** If the ribbon cable is not protruding, reach to the bottom of the *y* axis, stick in a finger, and push it out from within the *y* axis.

**Figure 10.** Connecting the injection unit ribbon cable

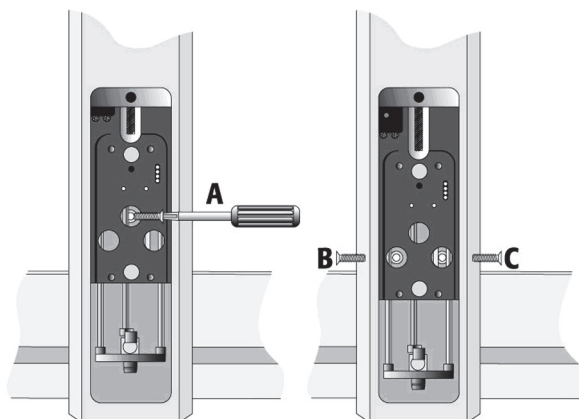


3. Hold the injection unit in place against the  $y$  axis. Make sure the two locating pins on the  $y$  axis fit into the two guide pin holes on the injection unit.
4. Place one of the screws onto the end of the supplied Torx driver. Slide the clear plastic cover on the injection unit all the way to the top. Locate the three large holes in the black anodized frame attached to the  $z$  axis inside the injection unit. Slide the frame upwards until the top hole is centered on the top threaded hole at the end of the  $y$  axis. Insert and securely tighten Torx screw A (Figure 11).

## 2 Installing the Accela Open Autosampler

Installing the Autosampler and DLW

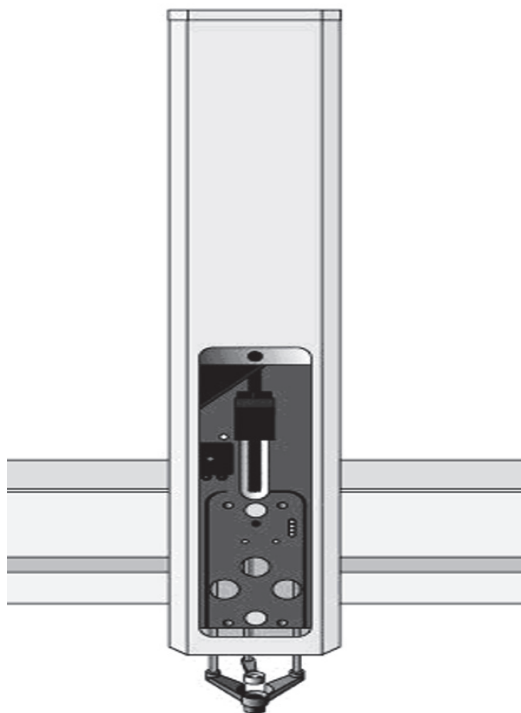
**Figure 11.** Inserting the injection unit mounting Torx screws



5. Install the two remaining Torx screws, B and C, in the left and right mounting holes, respectively. You might have to move the elastic cord slightly to the left to insert Torx screw C into the right hole.

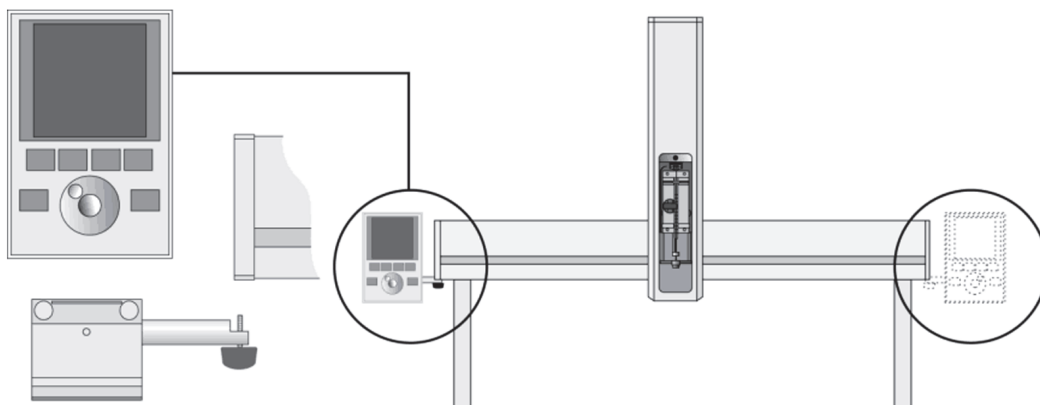
Figure 12 shows the completed installation.

**Figure 12.** Open Autosampler with injection unit



## Installing the Keypad Terminal

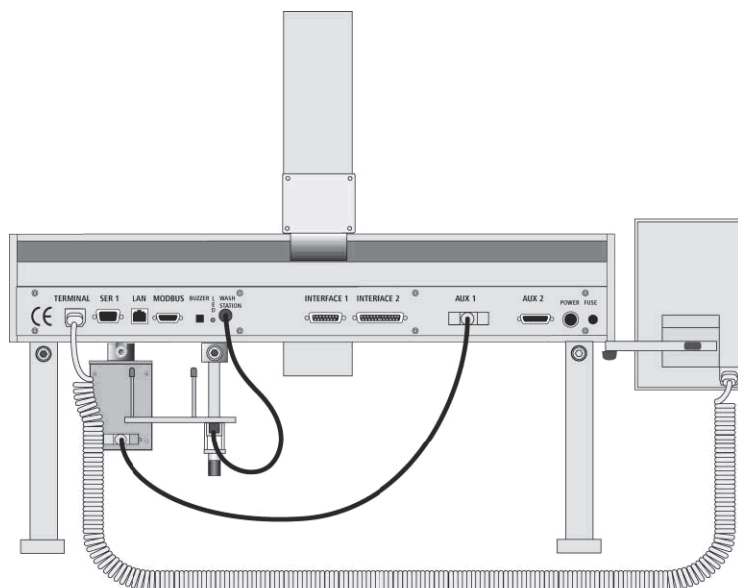
**Figure 13.** Installing the keypad terminal



### ❖ To install the keypad terminal

1. Install the safety shield on the left and right sides to the outside of the x axis. Use the provided, longer thumbscrew on the side where you plan to install the keypad.
2. Install the keypad mounting bracket on either the right or the left side of the x axis (Figure 13).
3. Connect one end of the white coiled cable to the keypad and the other end to the TERMINAL (SER3) interface jack on the back side of the x axis (Figure 14).

**Figure 14.** White coiled cable connection



4. Place the keypad terminal onto its mounting bracket.

## 2 Installing the Accela Open Autosampler

Installing the Autosampler and DLW

### Installing the Power Supply

#### ❖ To install the power supply

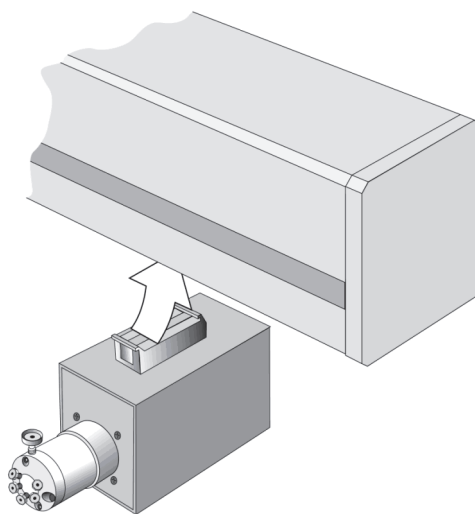
1. Locate the power supply, the DC power cable, and the AC power cable.
2. Set the power supply switch to OFF.
3. Connect one end of the DC power cable to the power supply and the other end to the POWER connector at the back of the *x* axis.
4. Connect the female end of the AC power cable to the power supply. Then connect the male end to an AC power outlet.

### Installing the LC Injection Valve

#### ❖ To install the LC injection valve

1. Locate the valve drive, which has one clamp that is identical to the other object clamps (Figure 15).

**Figure 15.** Installing the injection valve drive



2. Attach the valve drive to the *x* axis and tighten the mounting screw.

**Tip** To avoid delay volume between sample injection point and detection, place the injection valve near the detection device.

3. Connect the control cable from the valve drive to AUX1.

**Note** For information on installing another type or multiple valve drives, see “Electrical Connections for 2- or 4-Injection Valves” on page 37.

The injection valve and rotor are preinstalled on the valve drive.



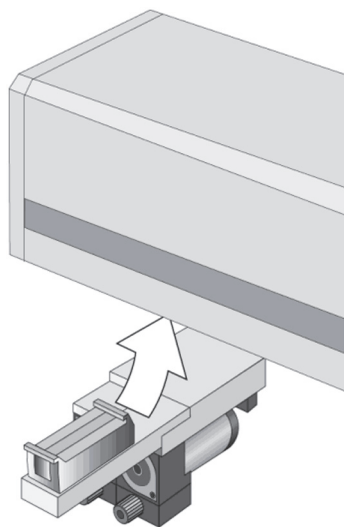
## Mounting the DLW Pump Holder

You must always mount the pump holder on the far left of the  $x$  axis. The solvent lines are connected to the holder. Moving across the pump holder with the  $z$  axis could kink the stabilizing wire. The best way to install the holder is by dismantling the drive from the Open Autosampler  $x$  axis.

### ❖ To mount the pump holder

1. Locate the pump holder and the two supplied Torx screws.
2. Attach the holder to the  $x$  axis as shown in [Figure 16](#).

**Figure 16.** Attaching pump holder bracket to the  $x$  axis



The solvent line assembly is not preinstalled at the factory. The correct mounting is shown in [Figure 17](#).

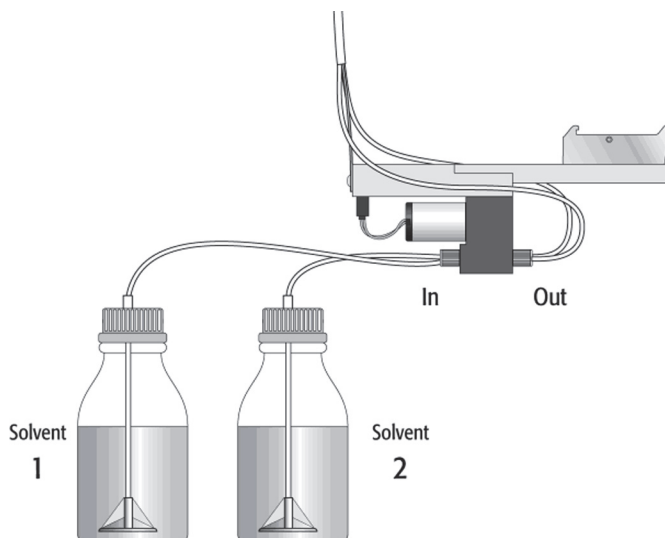
3. Using the end of the guide wire with the longer tubing ends, connect the guide wire with the screw (M4 x 8) and the corresponding serrated lock washer to the back of the module. Pay attention to the guide groove, which orients the wire ([Figure 17](#)).

**Tip** Always install the DLW tubing (from the kit) at the pump module first. Before you connect the other end of the tubing to the DLW Syringe assembly, make sure that the assembly is not installed in the injection unit.

## 2 Installing the Accela Open Autosampler

Installing the Autosampler and DLW

**Figure 17.** Installation of the solvent tube assembly to the pump holder

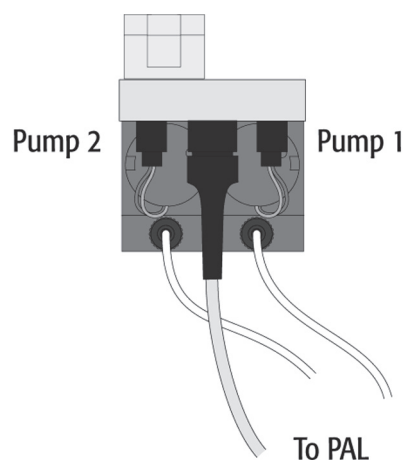


4. Connect the solvent tubing of the DLW tubing kit with the premounted nut to the pump module. Maintain the order: Solvent 1 connected to the left, and Solvent 2 to the right pump module as viewed from the front of the PAL System. This avoids any confusion when refilling or changing solvents.
5. Connect the solvent lines from the solvent reservoir bottle to the inlet connectors of the pump module.

The electrical connections from the pumps to the PCB sockets, mounted on the bracket, are done at the factory.

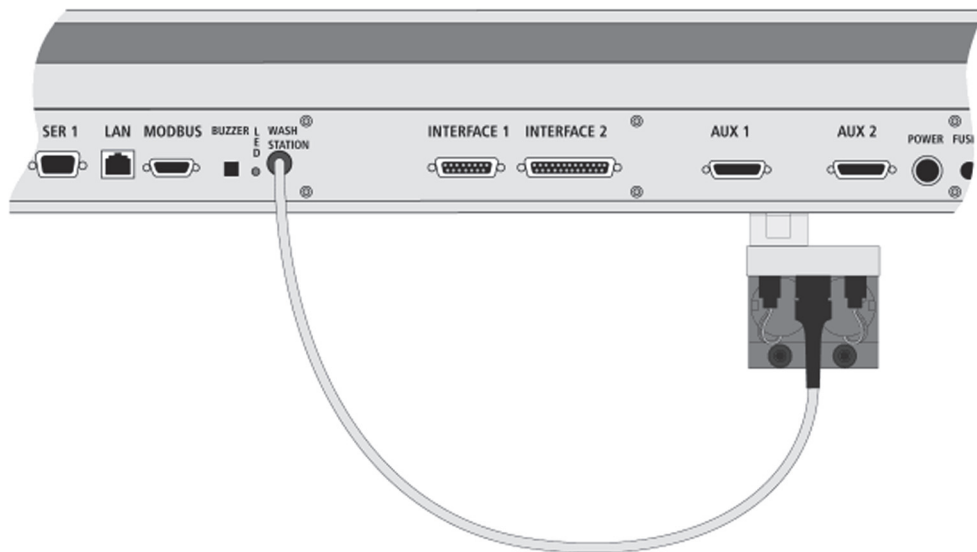
6. Insert the provided cable into the middle socket, between the other two sockets on the PCB. For details see [Figure 18](#).

**Figure 18.** Electrical connections at the pump holder bracket



The other end of the cable is connected to the wash station connector on the Control-xt board, as shown in [Figure 19](#).

**Figure 19.** Electrical connection from DLW pump holder to PAL-xt board



## Installing the DLW Option

This section provides instructions on installing the various parts of the DLW option.

- [Installing the DLW Wash Station](#)
- [Installing the DLW Syringe Holder Assembly](#)

### Installing the DLW Wash Station

You can connect the DLW Wash Station anywhere along the *x* axis. However, Thermo Fisher Scientific strongly recommends that you install it as close as possible to the injection valve. Make the travel path of the injection unit from the valve to the wash station as short as possible to keep the cycle time as brief as possible.

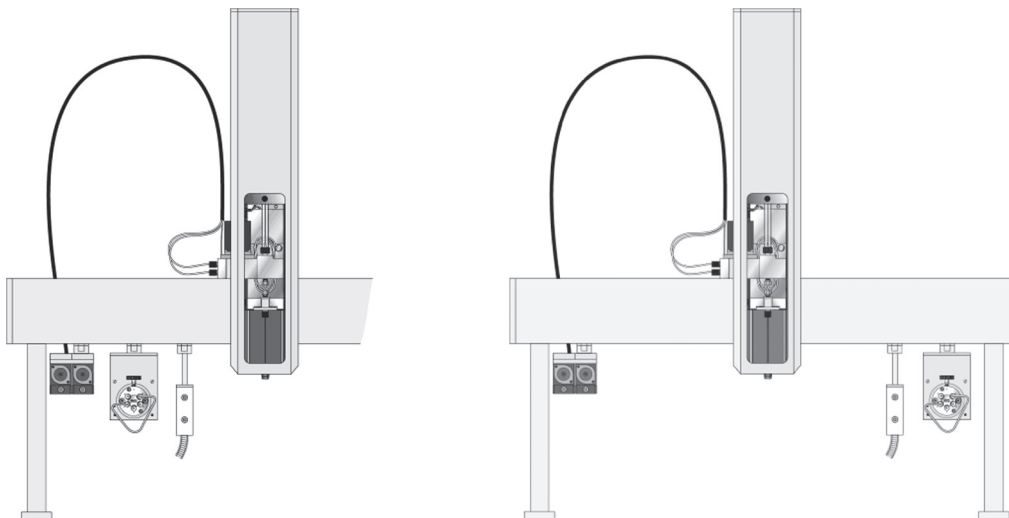
[Figure 20](#) shows two configurations for installing DLW modules, favoring a short connection from the injection valve to the other HPLC system components on the one hand, and always having the DLW Wash Station close to the injection valve on the other. Always attach the DLW Pump Module, with the connected DLW tubing from the kit, on the far left side of the *x* axis.

Such a configuration also has the advantage that you can connect the waste tube from the injection valve to the front waste inlet of the wash station module. This waste tube is part of the DLW option kit.

## 2 Installing the Accela Open Autosampler

Installing the Autosampler and DLW

**Figure 20.** Mounting the DLW option modules in combination with the injection valve



### ❖ To install the DLW Wash Station

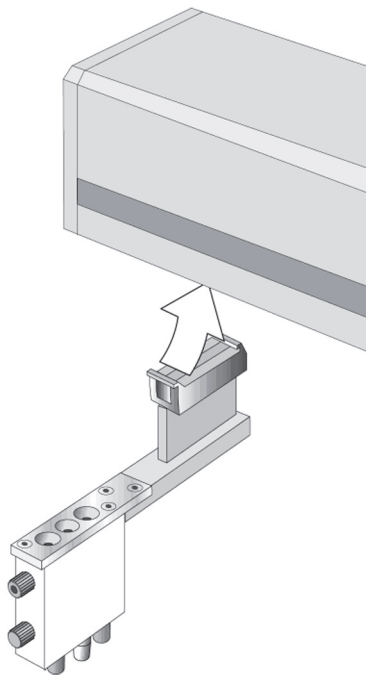
1. Locate the DLW Wash Station module.
2. Attach the holder to the  $x$  axis at the selected position as shown in [Figure 21](#).
3. Connect the waste from the injection valve to the DLW Wash Station.
4. Connect the waste tube to the wash station waste adapter.

Replace the dummy plug or plugs on the front if you are using them to connect the waste line coming from the injection valve. This simple connection of waste lines cleans up the system, avoiding several waste lines running to the waste container.

**Tip** Use the other two dummy plugs underneath the waste block for the DLW option.

You can connect the DLW tubing from the DLW pump outlet at these wash station inlet ports. The functionality of the fast wash station is then applied using this configuration.

**Figure 21.** Attaching the DLW Wash Station to the x axis



**Tip** The waste tube must be positioned below the injection valve. Make sure that the waste liquid flows into the waste container without restriction. The waste tube must always be above the liquid level in the waste container.

## Installing the DLW Syringe Holder Assembly

**Tip** Before you insert a syringe, first verify the position, Change Syr. Do not locate the Change Syringe position above an object where the needle could collide with a vial, for example.

Path: **Menu > Utilities > Syringe > function key F3 (Change Pos).**

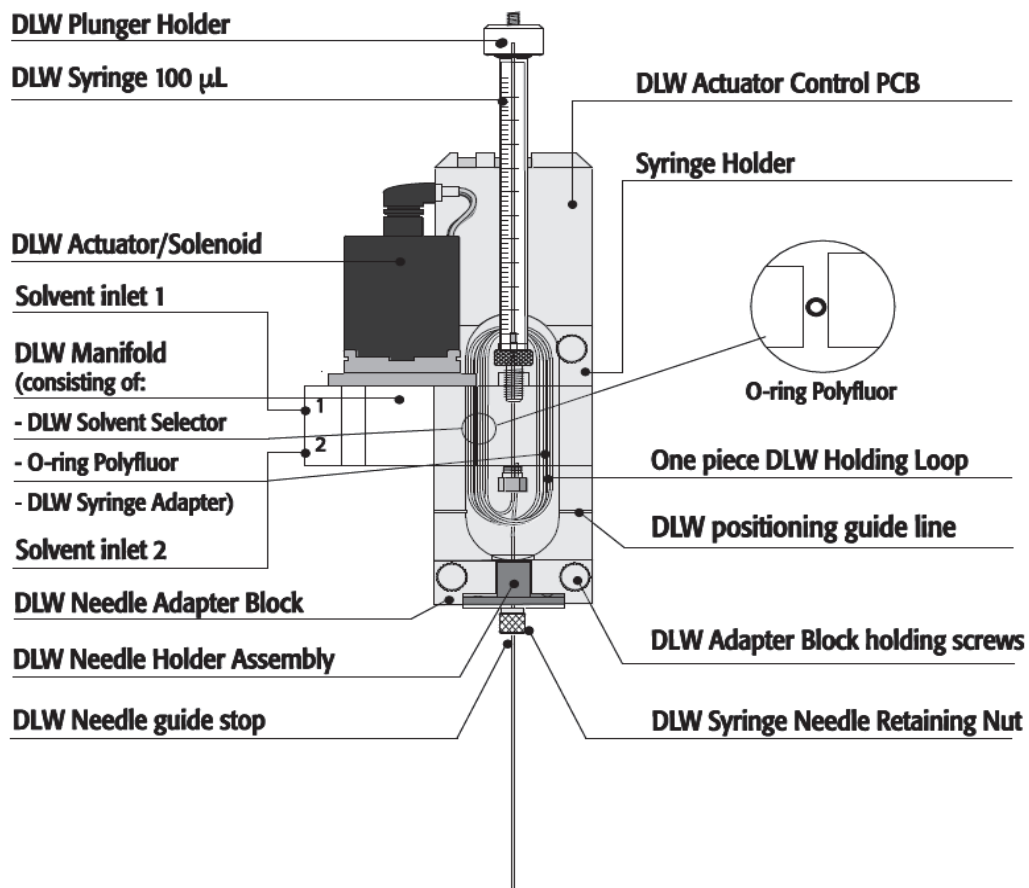
Move the PAL injection unit to a location where no collision can occur with the  $z$  axis by selecting the  $x$  and  $y$  axes appropriately. The value for the  $z$  axis is given as a default and a change of this position is not necessary in standard operation.

This precaution helps avoid needle damage during routine operation. Nevertheless, the description in the next section on how to install the DLW Syringe assembly recommends switching off the unit and inserting the assembly without using the Open Autosampler command, Change Syringe.

The DLW Syringe Holder differs from standard liquid syringe holders. The DLW manifold is attached to the syringe holder. The DLW Actuator/Solenoid and the Holding Loop connecting the syringe inlet (bottom) and Syringe Needle are attached to the DLW manifold. All of these parts are factory installed. See [Figure 22](#).

**Figure 22.** Front view of the DLW Syringe Holder assembly

### DLW Syringe Holder Assembly



To install the DLW syringe holder, follow these procedures:

- [Preparing and Installing the DLW Syringe](#)
- [Connecting the DLW Tubing to the Syringe Holder](#)
- [Inserting the DLW Syringe Holder Assembly into the Injection Unit](#)

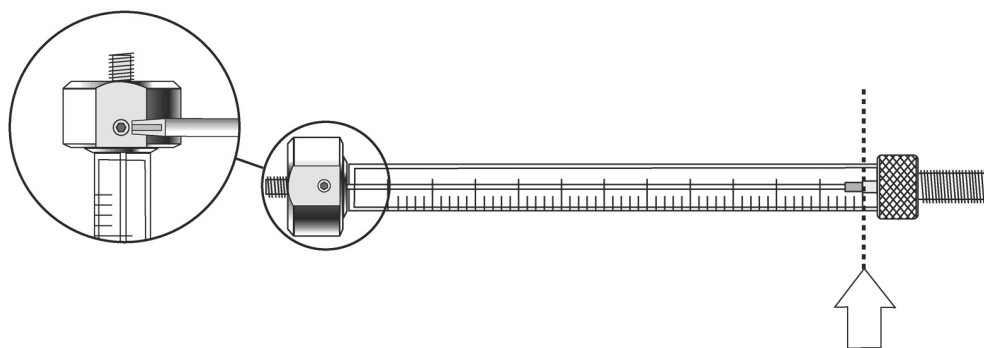
## Preparing and Installing the DLW Syringe

This section contains information on how to prepare and install the DLW Syringe.

### ❖ To prepare and install the DLW Syringe

1. Prepare the DLW Syringe by inserting the DLW Plunger Holder.
  - a. Move the plunger manually down to the stop position, and release a slight amount of pressure to the plunger tip by pulling a fraction of a millimeter backwards.
  - b. Install the DLW Plunger Holder and tighten the Allen screw (Allen key #6) firmly (Figure 23).

**Figure 23.** Inserting DLW Plunger Holder on DLW Syringe



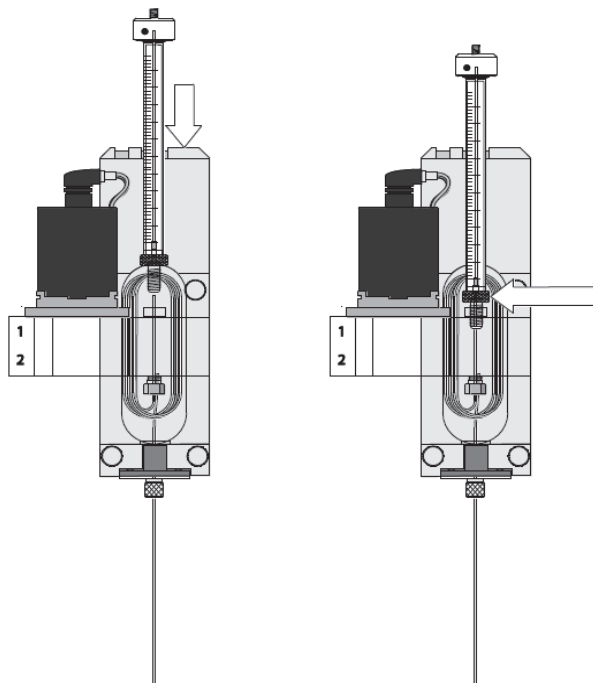
**Tip** Make sure the syringes are primed before beginning sample preparation. Prime all liquid syringes manually first before inserting them into the Open Autosampler system.

2. Screw the prepared DLW Syringe into the holder. Hold the syringe at the lower metal mount to tighten the syringe (Figure 24).

## 2 Installing the Accela Open Autosampler

Installing the Autosampler and DLW

**Figure 24.** Inserting DLW Syringe into holder



**CAUTION** Make sure to hold the syringe on the metal mount to tighten the syringe. Holding the glass barrel while tightening can damage the seal where the glass meets the metal.

**Note** Do not mount the syringe needle at this stage.

### Connecting the DLW Tubing to the Syringe Holder

**Note** Thermo Fisher Scientific recommends connecting the DLW tubing from the kit to the syringe holder while the assembly is not inserted in the  $x$  axis.

**Note** The wetted parts of the DLW Actuator/Solenoid are PEEK for the body and FFKM (Simriz) for the seal material.

PEEK tubing exhibits excellent chemical resistance to most of the chemicals used. However, the following solvents are not recommended for use with PEEK: DMSO, THF, methylene chloride (dichloromethane), nitric acid, or sulfuric acid. For more details, refer to the compatibility tables provided by the manufacturer of PEEK material or components.



❖ **To connect the DLW tubing to the syringe holder**

1. Connect the end of the sleeved tubing with the short tubes to the DLW Syringe Holder assembly. See “[Mounting the DLW Pump Holder](#)” on [page 23](#).
2. Use the groove in the block at the back of the DLW Manifold to orient the guide wire. To connect, use the provided screw (M3×5) and the corresponding serrated lock washer.
3. Test to see if the wire tension is enough to keep the tubing in an upright position but low enough to move the syringe holder manually along the *x* axis.

**IMPORTANT** The design of the DLW Manifold and the concept of washing by active pumping and active closing of the lines by the DLW Actuator theoretically do not require a predefined solvent line position.

A problem could occur if solvent is refilled or exchanged while the tubing still contains solvent from the previous setup. Always keep the lines at the same position, and prime the entire system with caution to prevent unnecessary confusion.

4. Connect solvent line 1 to the upper port and solvent line 2 to the lower port at the left side of the DLW Manifold.

Make sure the order of the tube connections, upper or lower, is consistent. In certain applications you cannot mix types of solvents; for example, biofluid sample solution should not come in contact with highly concentrated organic solvents.

**Inserting the DLW Syringe Holder Assembly into the Injection Unit**

**Tip** Insert the DLW Syringe Holder assembly when the Open Autosampler is powered off.

❖ **To insert the DLW Syringe Holder assembly into the injection unit**

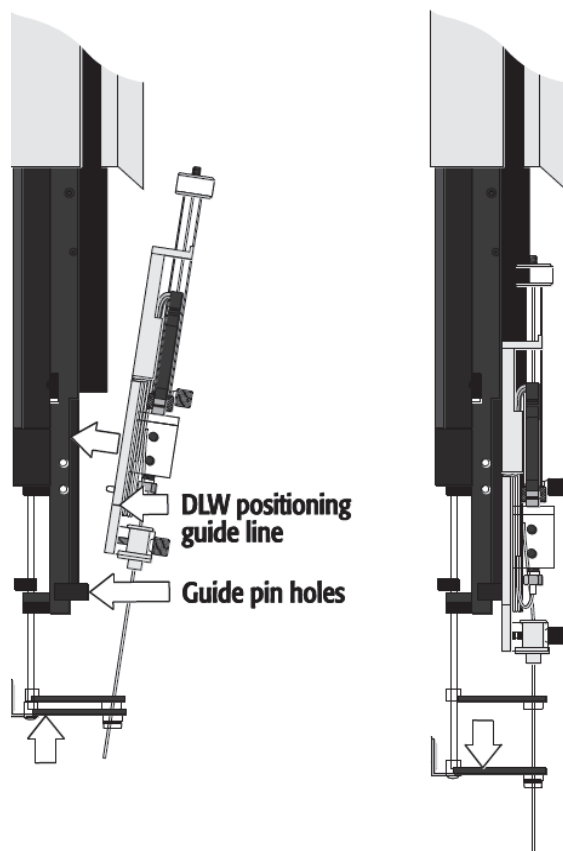
1. Manually move the PAL injection unit aside to allow free movement of the syringe slider. Lower the syringe slider ([Figure 25](#)) in order to gain access to install the DLW Syringe assembly.

Match the magnetic pins of the syringe holder using the counter positions at the syringe slider.

## 2 Installing the Accela Open Autosampler

Installing the Autosampler and DLW

**Figure 25.** Inserting the DLW Syringe Holder assembly into the injection unit



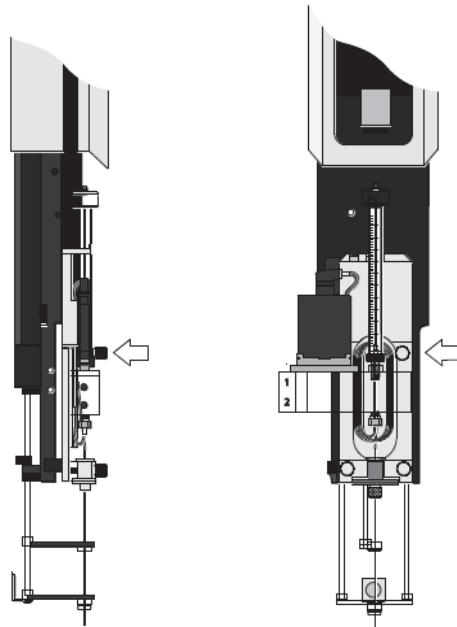
2. Press the syringe holder firmly against the z-axis slider to ensure that the holder engages.



**CAUTION** Do not press against the DLW Actuator/Solenoid. The mounting of the latter to the DLW Manifold is fragile and might break.

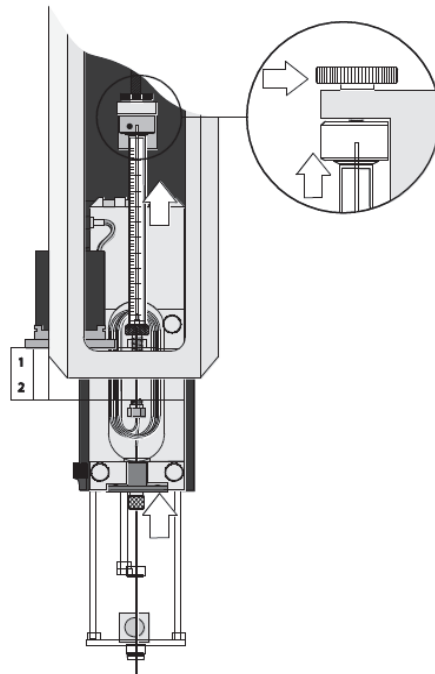
3. Tighten the knurled screw to fix the holder assembly to the syringe slider ([Figure 26](#)).

**Figure 26.** Fixing syringe holder to syringe slider



4. Move the plunger up (plunger holder) until the thread of the screw catches the thread of plunger bushing. Tighten the screw to fix the plunger holder ([Figure 27](#)).

**Figure 27.** Connecting the DLW Syringe plunger holder



## 2 Installing the Accela Open Autosampler

Installing the Autosampler and DLW

5. Tighten the holding screw to secure the syringe holder position.



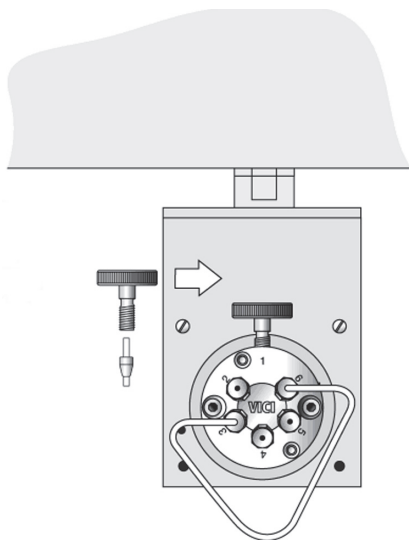
**CAUTION** Make sure the removable needle fits leak-tight to the syringe. A Teflon seal is provided. When installing the seal, ensure that the hole is open and that no fuzzy (lint) polymer material, which might clog the needle, is in the path. When slipping the Teflon seal over the needle, pay attention to the sharp edges of the needle that can cut into the polymer, which again might cause the needle to clog.

6. Move the lower needle guide carefully up and down to make sure that the needle tip does not catch on the guide.

### Installing the Valve Needle Guide Assembly

This section describes how to install the Valve Needle Guide assembly. Perform the installation as illustrated in [Figure 28](#).

**Figure 28.** Installing the standard valve needle guide



## Injection Valve Plumbing and Connection to DLW Wash Station

### ❖ To plumb the injection valve and connect it to DLW Wash Station

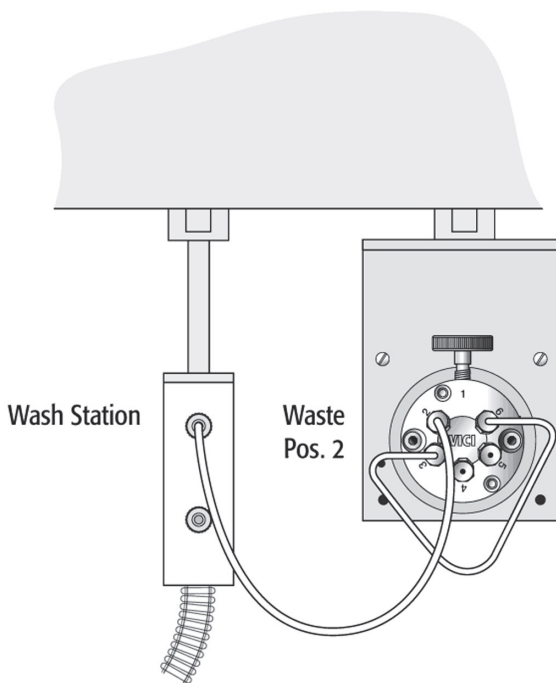
1. Connect the injection valve waste line from port 2 (for example, the standard Cheminert valve type), as shown in Figure 29, to the front waste port of the wash station.

**Tip** If you use two injection valves in the PAL configuration, connect the second waste line to the lower front waste port of the DLW Wash Station. If you use only one connector, apply a dummy plug to the second port on the wash station.

2. Place the waste tube from the DLW Wash Station into the waste container.

**IMPORTANT** You must position the waste container below the injection valve. Make sure that the waste liquid flows into the waste container without restriction.

**Figure 29.** Solvent line plumbing on Cheminert valve types



## 2 Installing the Accela Open Autosampler

Installing the Autosampler and DLW

### Electrical Connections

This section contains information on setting up the electrical connections of the Accela Open Autosampler.

#### Electrical Connections for Single Injection Valve Setup

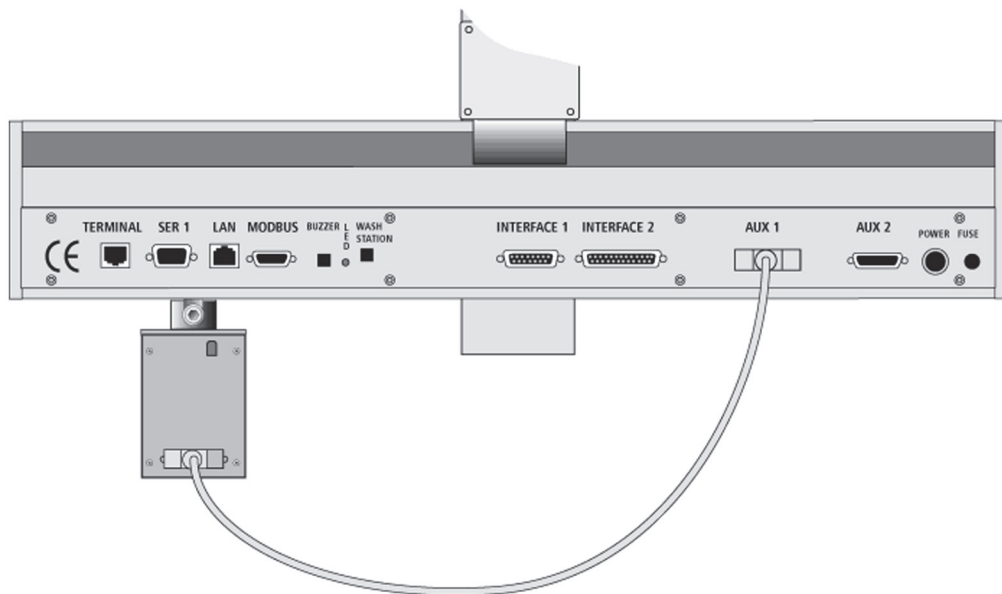


**CAUTION** Always power off the Open Autosampler before connecting or disconnecting the DLW cable or any other accessory cables.

The electrical connection of the DLW option is identical for the PAL HTS-xt and the HTC-xt models. Both models use the same board.

The diagram in [Figure 30](#) shows an installation using the valve drive connected to the AUX interface. If you use a serial or multiposition valve drive, see the diagrams for 2- or 4-injection valve combinations ([Figure 31](#), and [Figure 32](#) on [page 38](#)).

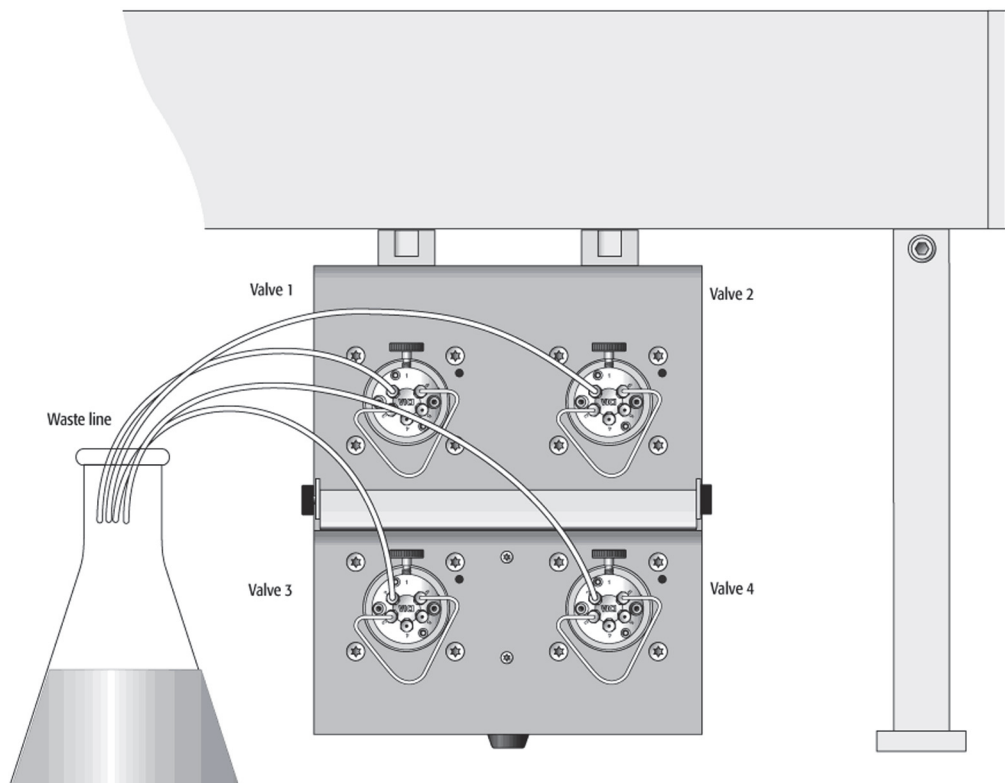
**Figure 30.** Electrical connection, DLW option on a PAL-xt



## DLW Option Installation in Combination with 2- or 4- Injection Valves

You can combine the DLW option with Open Autosampler systems using a setup of 2- or 4-injection valves (Figure 31). The concept of the DLW option is that the crucial parts of the wash station are attached to the injection unit.

**Figure 31.** DLW option in combination with 4-injection valves



The sample is taken up into the Holding Loop, and the syringe is actually the prepared reservoir for the following wash step after the loop filling and injection step.

Following this concept allows for more than one injection valve in the Open Autosampler configuration. Neither a special setup nor additional modules are necessary to inject into and clean one or more valves connected to the Open Autosampler system. As mentioned, the crucial DLW parts are attached to the injection unit at any available position.

## Electrical Connections for 2- or 4-Injection Valves

Connect the valve drives as shown in Figure 32.

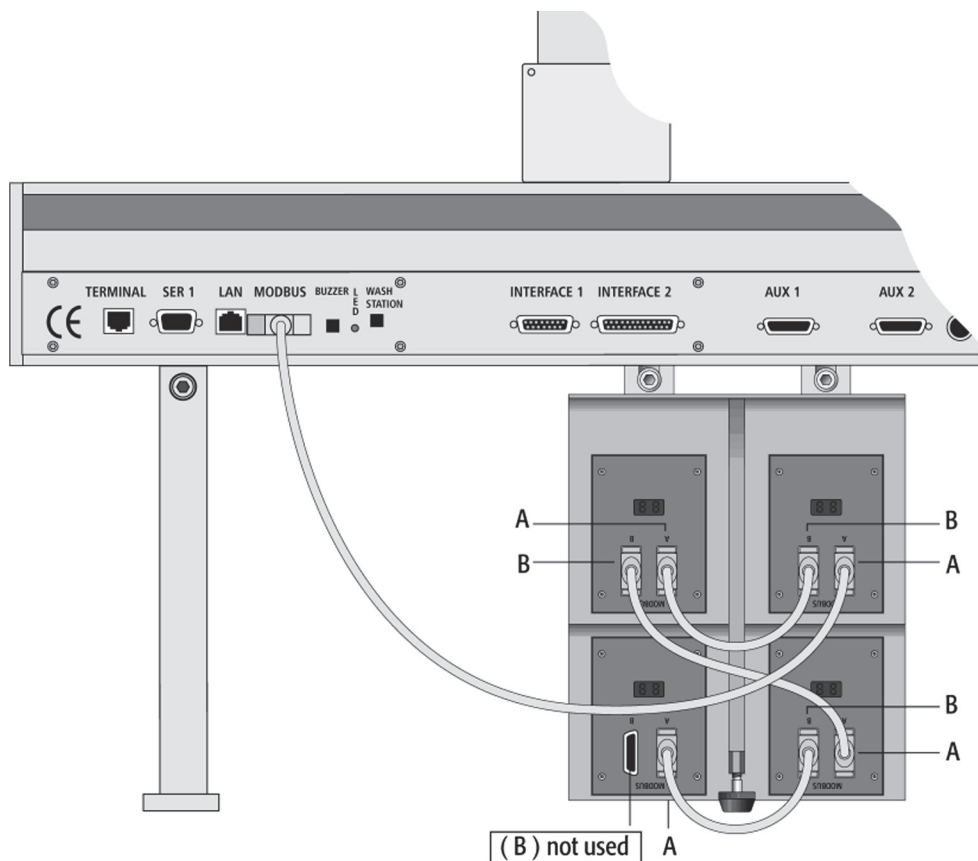
Because you can only use the DLW option in combination with the PAL-xt systems, you must connect the first cable from the PAL Interface MODBUS to the serial valve drive connector A. Each drive is connected with daisy-chain cabling.

The last serial drive connector B remains open or unused. The interface MODBUS provides the electrical current and interface logic.

## 2 Installing the Accela Open Autosampler

Installing the Autosampler and DLW

**Figure 32.** Electrical connections to 4-injection valves, daisy-chaining with PAL-xt system



### Installing the Power Supply for the Cooled Stack

The power supply for the cooled stack has a built-in control unit with display (P/N CH-952995).

#### ❖ To install the power supply

1. Locate the power supply and the AC power cable.
2. Set the power supply switch to OFF.
3. Connect the open end of the DC power cable from the stack/tray cooler to the power supply connector labeled, Peltier Thermostat.
4. Connect the female end of the AC power cable to the power supply, and then connect the male end to an AC power outlet.
5. Set the switch on the power supply to ON.
6. Observe the display to ensure that it reads the preset temperature of +10 °C.



## System Synchronization Connections

The system interconnect cables and adapter cables that synchronize the run signals for an LC or LC/MS system with an Accela Open Autosampler depend on the Accela pump model, the mass spectrometer model, and whether the LC system includes an Accela detector.

Table 5 lists the system interconnect and adapter cables that Thermo Fisher Scientific supplies with the Accela Open Autosampler or by special order.

**Table 5.** Accela Open Autosampler contact closure and adapter cables

Cable description	Part number	Use
Accela Open Autosampler system interconnect cable  (supplied with autosampler)	60157-63024	Required for all hardware configurations
Accela detector and MS detector adapter cable  (supplied with autosampler)	60157-63026	Required for hardware configurations that include one or more of the following: <ul style="list-style-type: none"> <li>• Accela detector</li> <li>• TSQ or Exactive mass spectrometer</li> </ul>
Accela Pump adapter cable  (available by special order)	60157-63022	Required for the Accela Pump

**Note** Thermo Fisher Scientific has discontinued the Accela Pump.

For instructions on how to connect the required cables listed in Table 5, refer to the *Accela Open Autosampler User Guide*.

# Accela Open Autosampler Object Positions

This section contains information about Accela Open Autosampler object positions.

## Defining Object Positions

**Note** Remove the syringe adapter from the injection unit before performing the following steps.

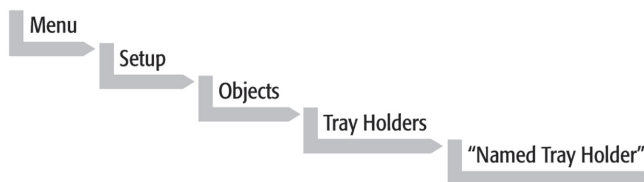
The objective is to define the reference positions for all Accela Open Autosampler objects. Make sure to properly mount the tray holders, valve drives, and wash station to the Accela Open Autosampler  $x$  axis. The following describes how to teach the reference position for a tray holder. The described procedure is common to all Accela Open Autosampler objects.

### ❖ To teach the reference position for a tray holder

1. Switch on the Accela Open Autosampler power supply.
2. Observe the keypad display.

The model name, Accela Open Autosampler, appears along with the firmware version number. The Job Queue menu screen then appears.

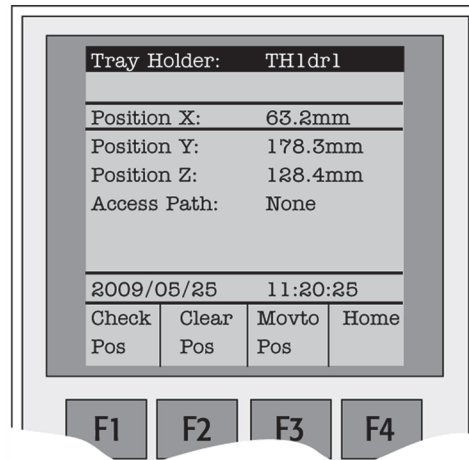
3. Complete the following sequence (common to all objects):



where the Named Tray Holder represents a predefined tray holder (for example, THldr1 or Stack1).

After selecting Named Tray Holder, the  $x$ ,  $y$ ,  $z$  positions associated with the object are displayed (Figure 33).

**Figure 33.** Menu screen object Tray Holder



4. Select **Position X** and press ENTER.

The injection unit moves to the previously defined *x*-axis position.

5. Rotate the outer knob to adjust the *x*-axis position to the tray holder reference position.
6. Press the inner knob to enter the Position X value.
7. Repeat [step 4](#) through [6](#) for Position Y and Position Z.

If the F3 button (Move to Zero) is activated, the injection unit moves to the HOME position.

8. Verify the defined *x*, *y*, *z* positions by pressing F1: Check Pos.

The release of PAL firmware level 4.1.2 makes possible the correction of all three axes, *x*, *y*, and *z*, in the firmware class tray. You can correct the inclination for a tilted tray in the direction of the row, the column, or both. Teaching is possible in the Utilities section by using the path: **Menu > Utilities > Trays**.

## Description of Object Positions

This section contains a description of the object positions.

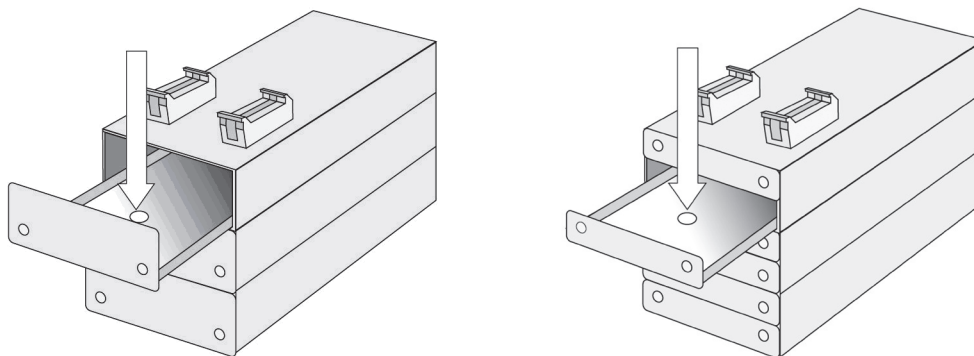
### Tray Holders Stack or Stack Cooler with 3 or 6 Drawers

The reference position for a stack or stack cooler is identical. The reference position in the case of a 3-drawer stack is a hole in the upper drawer and in a 6-drawer stack, in the second drawer from the top (same height level), as shown in [Figure 34](#).

## 2 Installing the Accela Open Autosampler

### Accela Open Autosampler Object Positions

**Figure 34.** Stack reference position



For the object stack/tray cooler, you must properly position the  $x$ ,  $y$ ,  $z$  axes in the Open Autosampler system.

After  $x$ -,  $y$ -,  $z$ -axis positioning of stack cooler model 6DW or 12MT, add the following steps to verify the correct installation of the module.

❖ **To verify the correct installation of the module**

1. Close the drawer manually.

Choose **Menu > Setup > Objects > TrayHolders > *corresponding device***.

Press function key F1, Check Pos. The injection unit moves to the predefined position.

2. Open the bottom drawer manually.
3. Remember the  $z$  value. Turn the  $z$  axis all the way down to check that the black needle guide of the injection unit fits neatly into the position hole of the drawer.
4. If the position is proper, turn the injection unit slider back to the original  $z$  value. Then close the drawer.
5. If the position is off by one or two millimeters, compensate for any discrepancy in the  $x$ ,  $y$  position seen between the upper and lower drawer.
6. If the position is off by more than one or two millimeters, check the bracket attachment of the module to the  $x$  axis carefully. If in doubt, loosen the claws again and reposition the unit on the  $x$  axis.

If this action is not successful, check the  $x$ - to  $y$ -axis angle and the  $y$ - to  $z$ -axis angle. If one of these angles is not  $90^\circ$ , make corrections.

Ensure also that the  $y$  axis runs parallel to the open drawer. If this is not the case, the bracket alignment is probably out of square.

**Note** The  $y$  axis tolerance for the teaching point is limited to 129.4 mm (5.09 in.).

When defining the position of the stack cooler, make sure you stay within this tolerance with the reference point at the top drawer.



**CAUTION** Do not use vials without a sealing cap or a microtiter- or deepwell-plate without a plate seal. Vapor phase from organic solvents can be hazardous and flammable. Acidic vapor phase can cause corrosion to critical mechanical parts.

## Trays

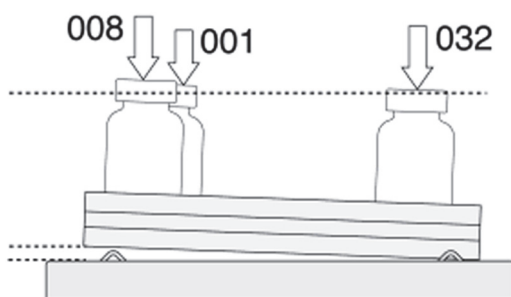
Generally, you do not need to teach a tray position. The tray holder has a teaching position, as described earlier. The firmware assigns a tray to the tray holder. The numbers (values) for the relevant items are stored in the firmware object tray. To complete the combination of a tray holder and a tray, you must assign the corresponding tray type to the tray.

The tray type contains the geometric data for the tray and information, such as how many samples are in a row and a column in the tray. A tray type defines the pattern and sampling sequence of the sample location within a tray.

Again, the tray is physically placed in the tray holder, the firmware assigns the tray to the tray holder and the tray type to the tray. This is routine procedure as long as no special circumstances intervene, such as customized trays or tray types.

PAL Firmware level 4.1.2 or later now makes a correction possible if the tray surface is not exactly horizontal and planar but is inclined in one or more axes. The ability to make the correction is standard at level 4.1.x. A dialog box opens the moment the tray has been checked out at the corners to verify vial positions and heights. The following sections explain how to correct a possible inclination ([Figure 35](#)).

**Figure 35.** Demonstrating a possible inclination of the tray in  $x$ ,  $y$ ,  $z$  axes

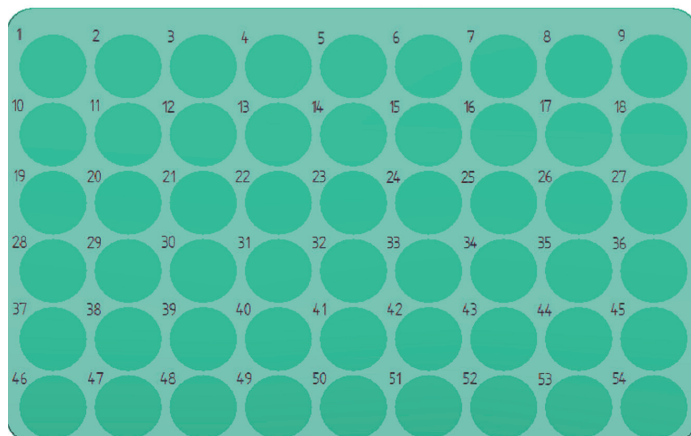


### Definition of a Tray Row and Column

The Accela Open Autosampler defines rows and columns by the order in which samples are treated. A row is not associated with an  $x$  or  $y$  axis. [Figure 36](#) provides a visual definition with tray type VT54.

For additional information about sample trays, refer to the *Accela Open Autosampler User Guide*.

Figure 36. Tray type VT54

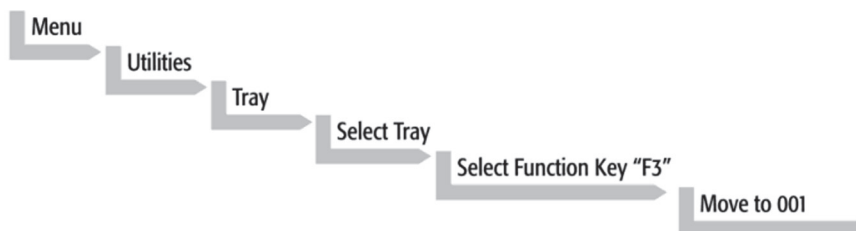


### Correction for Inclination in X, Y, or Z Axes

The following describes the correction using the path to the Utilities section.

#### ❖ To correct for inclination in x, y, or z axes

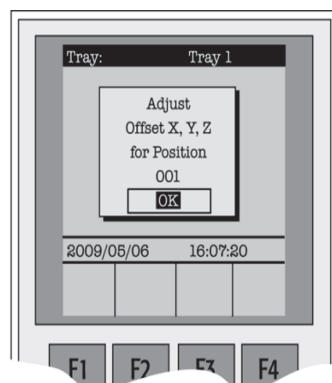
1. After selecting the Utilities function, Tray, verify whether the tray type assigned corresponds to the tray placed into the tray holder. If so, check the vial positions at the three corners of the tray. The F3 function key activates this.



A message box opens, displaying a note: Adjust offset X-, Y-, Z for Position 001.

2. Select **OK** to adjust all three axes perfectly by the top of the vial. This step is only valid for position number 001. See [Figure 37](#).

Figure 37. Message to adjust offset x, y, z for position 001

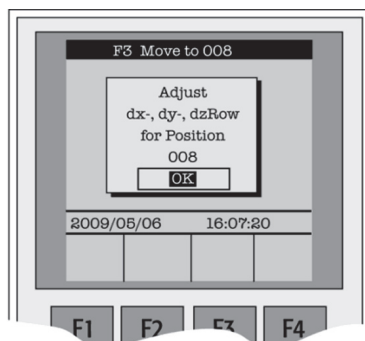


3. To check the next corner of the tray, activate the F3 function key again, for example, tray type VT54, position 009.

The message that appears prompts you to correct a possible inclination of the row.

4. Select **OK** to adjust the three axes perfectly to match the vial top or, in case of a well-plate, the top of the plate. See [Figure 38](#).

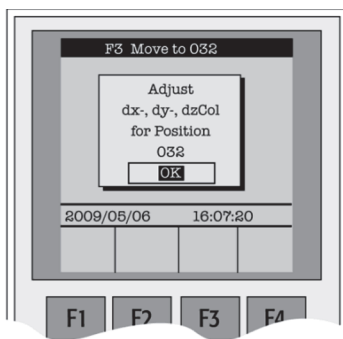
**Figure 38.** Message to adjust row inclination at the second tray corner position



5. To move to the next corner of the tray, activate the F3 function key for position 54 in the example of tray type VT54.

You can correct a possible inclination of the tray column in the same manner as you did with the first two corners. See [Figure 39](#).

**Figure 39.** Message to adjust column inclination at the third tray corner position



The three corner points are now adjusted for a possible inclination of the tray in any axis. Without further teaching, the Accela Open Autosampler interpolates a possible deviation from an ideal axis position for the other vials, caused by variance from the horizontal.

**Note** For routine work using a standard vial, you do not need to compensate for a possible inclination as long as the vial detection is within the Z-Tolerance range.

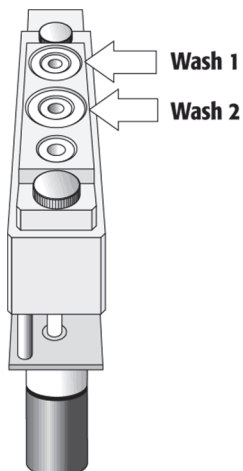
## 2 Installing the Accela Open Autosampler

Accela Open Autosampler Object Positions

### Wash Station: Wash1/Wash2 Reference Point

Figure 40 shows the Wash Station Wash1/Wash2 reference point.

**Figure 40.** Wash station Wash1/Wash2 reference point

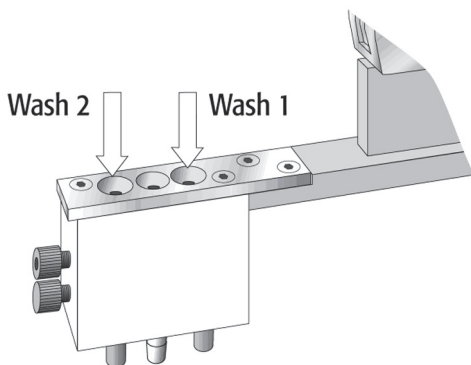


For a fast wash station, the reference positions are the two holes in the wash vial caps (Figure 40). Be sure to center the lower needle guide in these holes with the bottom of the lower needle guide lightly touching the surface of the wash station assembly.

For the DLW Wash Station option, the reference positions for Wash1 and Wash2 are the two holes on top of the wash station assembly. See Figure 41.

Center the lower needle guide in these holes with the bottom of the lower needle guide lightly touching the surface of the wash station assembly.

**Figure 41.** DLW Wash Station Wash1/Wash2 reference point





❖ **To make the reference position of Wash Station Wash1/Wash2**

1. Press HOME on the handheld controller and select the sequence:

**Menu > Setup > Objects > Wash Stations > Wash1**

2. Enter **Wash1** and adjust the  $x$ ,  $y$ ,  $z$  positions until the lower needle guide of the injection unit centers on top of the wash station assembly. Fine-tune the  $z$  position until the bottom of the lower needle guide is flush with the top surface.

3. Repeat steps 1 and 2 for Wash2 by using this path:

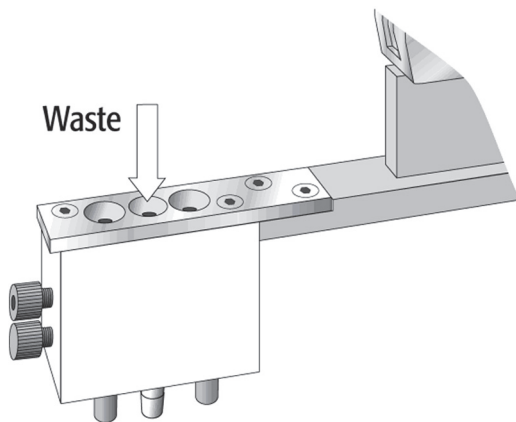
**Menu > Setup > Objects > Wash Stations > Wash2**

**Waste Reference Point, Injectors (Waste)**

The waste position represents an injector within the Accela Open Autosampler software. It is defined in the object class, Injectors.

The reference position for the waste port—a hole (slightly larger than the needle guide)—has been placed in the middle position of Wash1 and Wash2 (Figure 42). Center the lower needle guide in this hole with the bottom of the lower needle guide lightly touching the surface of the wash station assembly.

**Figure 42.** Waste reference position



❖ **To make the Waste reference position**

1. On the handheld controller, select this sequence:

**Menu > Setup > Objects > Injectors > Waste 1**

2. Select **Waste** and adjust the  $x$ ,  $y$ ,  $z$  positions until the lower needle guide of the injection unit centers on top of the wash station assembly. Fine-tune the  $z$  position until the bottom of the lower needle guide is flush with the top surface.

3. Press F1 to verify the position and F4 for Home.

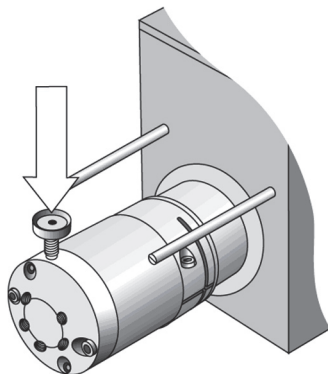
## 2 Installing the Accela Open Autosampler

Accela Open Autosampler Object Positions

### Injection Valve Reference Points

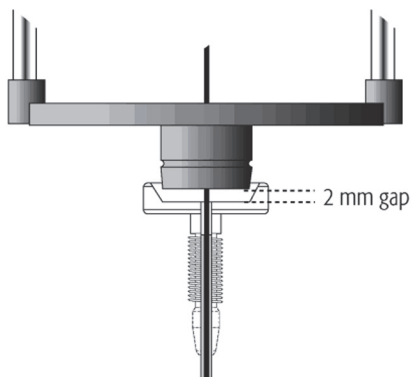
For an LC Valve the reference position is the valve needle guide fitting mounted on the top valve port (Figure 43).

**Figure 43.** LC Valve reference point



The lower needle guide of the injection unit is centered in the valve needle guide fitting. Adjust the z position so that the bottom of the lower needle guide just touches the surface of the valve needle guide fitting. Then reduce the value by 2.0 mm. See Figure 44.

**Figure 44.** Positioning z-axis needle guide on valve needle guide



### Defining the Valve Penetration Value

Before performing the following steps, make sure object positions X, Y, and Z for LC Vlv1 are properly defined and the complete DLW syringe assembly is inserted into the injection unit.

#### ❖ To define the valve needle penetration depth

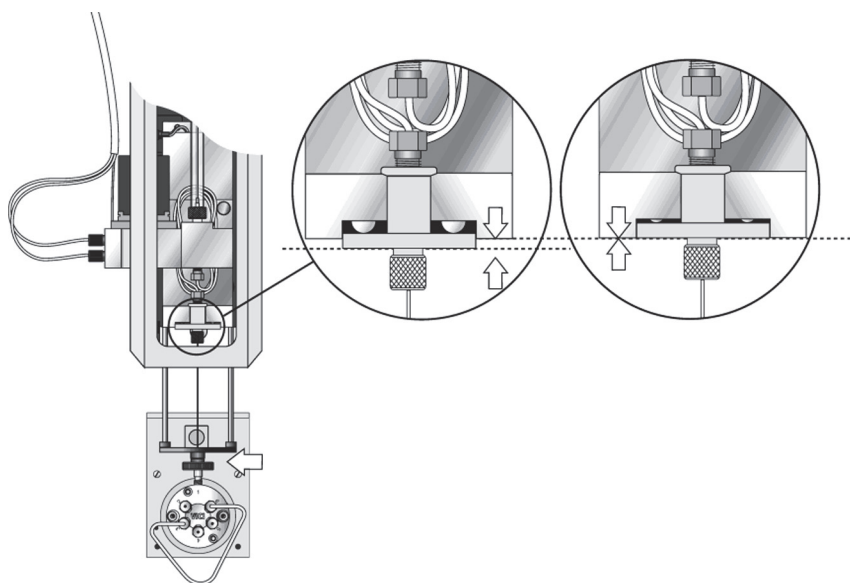
1. On the handheld controller select this sequence:

**Menu > Setup > Objects > Injectors > LC Vlv1**

2. Press F1, Check Pos. The injection unit moves to the previously defined LC Vlv1 position.
3. Select the item **Needle Penetr** with the cursor bars by pressing ENTER.

4. Slowly rotate the outer knob to adjust the needle penetration depth. The syringe moves down stepwise into the injection port.
5. When the syringe needle tip enters the valve needle guide, slow down the Z-movement again. Always observe the syringe needle during this adjustment step.
6. Move down the syringe stepwise until the bottom of the cross bar of the DLW Needle holder assembly is flush with the lower line of the DLW needle adapter block. See [Figure 45](#).
7. Press ENTER to save the needle penetration depth value.
8. Verify the defined needle penetration depth value by repeating steps 3 through 6.

**Figure 45.** Teaching the needle penetration



9. If you hear a click and the ball-spring loaded needle holder is completely compressed, stop the Z-down movement immediately.
10. Rotate the outer knob stepwise in the opposite direction until the gap of the spring loaded needle holder is approximately 1 mm, or the bottom of the holder assembly is flush with the lower surface of the needle adapter block, as in [Figure 45](#). Repeat and verify the needle penetration test.

[Figure 46](#) shows an example of the correct and incorrect positioning of the syringe needle tip in the valve inlet port.

**Figure 46.** Valve needle penetration depth



The ball-spring loaded needle holder presses the square cut needle tip firmly against the bottom of the valve body. This ensures a constant seal during the injection process.

## Injection Valve

This section contains information about the injection valve.

### Valve Drives and Valves – General Remarks

An Accela Open Autosampler can be equipped with one LC injection valve connected and controlled through the auxiliary interface (AUX). If you must configure more than one valve, or if an auxiliary interface is occupied by another module, such as a dilutor, then the autosampler requires multiposition or serial valve drives. The serial valve drives are connected at MODBUS interface.

The Accela Open Autosampler is equipped with the control board APR Control-xt, which provides two auxiliary interfaces, AUX1 and AUX2. You have the choice to connect one more valve drive.

### Injection Valve Flow Path

The six-port valves are connected to the detection system according to the flow paths as shown in [Figure 47](#).

#### Standard LC Injection Valve, Cheminert 6-Port

The standard LC injection valve is a six-port, two-way switching valve with a flat rotary plate and has a special needle guide ([Figure 47](#)).

**Figure 47.** Flow path of standard LC injection valve (Cheminert type)

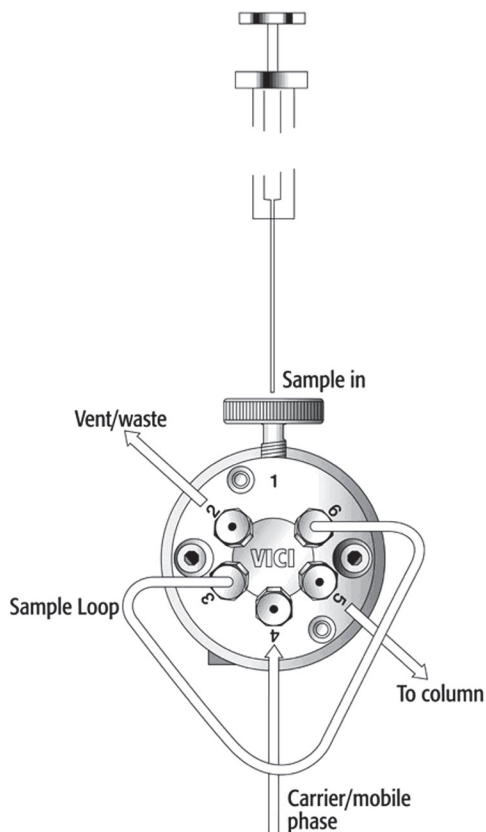


Table 6 lists the specifications for a standard LC injection valve.

**Table 6.** Standard LC injection valve specifications

Component	Specification
Valco/VICI Cheminert	Flat plate rotary valve
Stator material	Stainless steel 316
Rotor material	Valcon H (inert mix-polymer)
Port size	1/16 inch
Fittings	Valco 1/16 inch ferrules and nuts
Actuator	PAL valve drive (micro-electric actuator)
Loop size	20 $\mu$ L (shipped), other loop sizes available
Bore size	0.25 mm
Port to port volume	320 nL for vertical port, 75 nL for standard ports, 75 nL for engraving

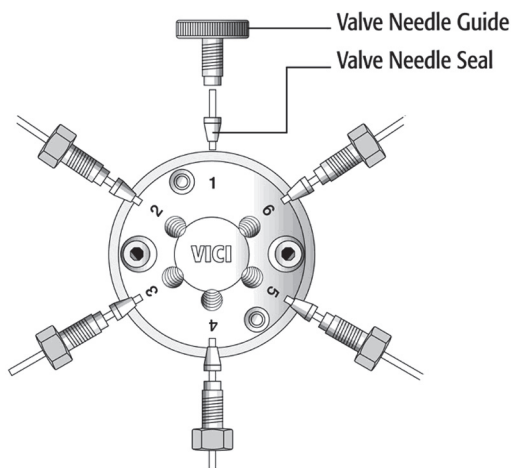
**Note** Sample loops of the different valve types are not interchangeable.

## Valve Needle Guide and Needle Seal

All valves are equipped with a special valve needle guide fitting. This fitting has a wide diameter to connect with the syringe needle guide on the injection unit. The valve needle guide also has a countersunk hole to guide the syringe needle into the injection port.

Figure 48 shows an example of the setup with the standard valve, Cheminert type.

**Figure 48.** Valve needle guide and valve needle seal



The valve needle guide holds the needle seal. The needle seal, which is a short length of FEP tubing, forms the seal around the syringe needle. A stainless steel ferrule is tightened around the FEP sleeve to ensure a leak-proof fit.

### ❖ To ensure reproducible sample injection and minimize carryover

1. Properly install the valve needle guide and the valve needle seal.
2. Change the valve needle seal at regular intervals.
3. Set the needle penetration depth accurately (“[Injection Valve Reference Points](#)” on [page 48](#)).

To specify a maximum number of penetrations into the needle seal is not practical. You might want to check the tightness of the seal daily or weekly when first using the Accela Open Autosampler to gain experience with the specific use in the laboratory related to the application.

### ❖ To test the tightness of the seal

1. Take the needle seal out of the valve port.
2. Manually insert a syringe needle gauge 22 (or gauge 22S) from the top, down into the tube, and past the sealing point of the ferrule.
3. If you observe a restriction, continue to use the needle seal. If not, replace the needle seal.

## Interval to Replace Needle Seal

The interval to replace the needle seal depends on the number of penetrations. Consider these two factors when deciding to replace the needle seal:

- Leak, seal not reliable

Remove the needle seal from the inlet port and insert a syringe manually with the corresponding needle gauge (standard Gauge 22 OD 18/25 mm). Move the needle below the seal point of the ferrule. If you do not observe resistance at this point, replace the needle seal.

- Carryover

If you suspect any contamination of the needle seal with a compound, replace the needle seal.

Contamination can occur if the sample solution is accidentally dispensed too fast into the valve system. The restriction of valve and loop might cause backpressure, forcing the sample solution into the inlet port of the valve.

**Note** PAL Firmware (level 2.4.x or later) provides a counter for the number of injections performed using the PAL System. This counter monitors the valve switches and not the number of penetrations into the injection port; for example, cleaning the valve requires an extra penetration. You might want to monitor this counter in conjunction with the test for the tightness of the needle seal as described above.

After the autosampler runs routinely for a few weeks, the warning limit for the inject counter gradually becomes evident, making it an invaluable tool in routine work.

Choose **Menu > Info > Maintenance > Inject Limit**. (Inject counts reflect the actual number of injections.)

Table 7 lists the part numbers for replacement needle seals.

**Table 7.** Part numbers for replacement needle seals

Part	Description	Part number
Standard needle seal, pkg of 10	FEP Tubing transparent, ferrule SST, needle gauge 22 (18/25 mm OD), for Valco/VICI SST valves, Cheminert and W-type	CH-952451
Needle seal, pkg of 10	FEP Tubing transparent, ferrule PEEK., needle gauge 22 (18/25 mm OD), for Valco/VICI PEAK valves, Cheminert-type	00950-01-00331





# Operating the Accela Open Autosampler

This chapter describes how to operate the Accela Open Autosampler including how to use the control terminal, methods, job queue, utility functions, and logfile; and how to set up functions.

## Contents

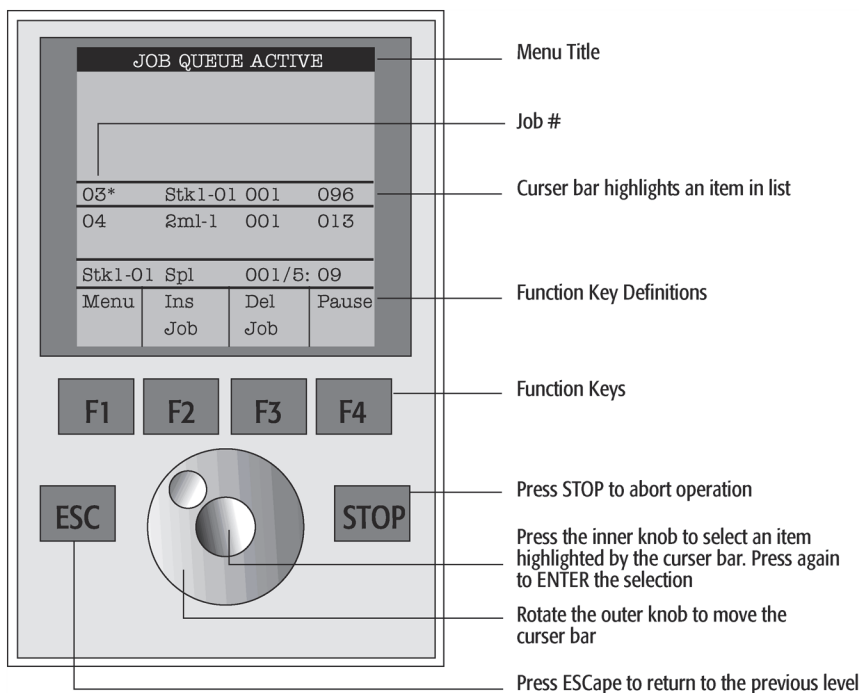
- [Control Terminal](#)
- [Methods](#)
- [Job and Job Queue](#)
- [Utility Functions](#)
- [Logfile](#)
- [Info Functions](#)
- [Setup Functions](#)
- [Communicating with the Computer](#)

## Control Terminal

The following procedures present the key steps required to set up and process multiple groups of samples with the Accela Open Autosampler. These procedures provide an overview for new users and a reminder for infrequent users. Make sure that the Accela Open Autosampler and all accessories are installed with objects defined correctly. In addition, ensure installation of a syringe of the specific type called for by a particular method.

[Figure 49](#) illustrates the Open Autosampler control terminal and the conventions used to enter, edit, and view information.

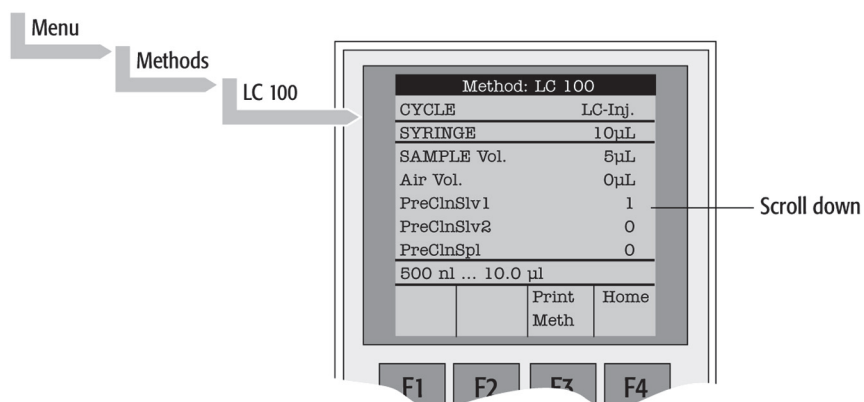
**Figure 49.** Accela Open Autosampler control terminal and conventions



## Menu Screens

The control terminal displays different menu screens, depending on the Accela Open Autosampler operating status and the particular function that you are accessing. All menu screens have the same basic format. The menu title displays at the top of the screen. A list of items displays below the title. The date and time, or status, show in the highlighted area above the function key labels on the bottom of the screen (Figure 50).

**Figure 50.** Accessing a method screen



## Function Keys

Functions (keys F1, F2, F3, and F4) correspond to the options for a particular menu, directly above each function key label.

Pressing the function key labeled Home always returns to the Job Queue menu.

## ESC and STOP Keys

Press ESC to return to the previous menu. Press STOP to cancel the current cycle, job, or job queue.

## Scroll Knob and Enter Button

Rotate the outer knob to scroll through items in a menu list. To select a highlighted item, press the central knob (the Enter button). Then use the outer knob to scroll through available options for that item or to change a numeric value. Press the inner knob again to enter the displayed option. You can also use the inner knob for other operations that require an Enter operation to continue or complete an operation.

## Methods

This section contains information about how to create, edit, view, and delete methods.

### Creating Methods

You can define methods and assign them names up to eight characters in length. You can create, copy, edit, and view methods from the Methods menu. Although you can view methods from the Job Queue menus, you cannot edit them there.

You create methods by either copying an existing method or creating a new method.

#### ❖ To copy an existing method

1. Choose **Menu > Methods > *method name* > Copy Method**.
2. At the prompt, enter a name for the new method.
3. Use the scroll knob and the left-right arrow function keys (F2 and F3) to select among alphanumeric characters and spaces.
4. Press the ENTER key to accept the name.

#### ❖ To create a new method

1. Choose **Menu > Methods > Select the Insertion Point > Instrument Method**.
2. Enter a new method name.

After you create a copy of the method, the method parameters appear and can now be edited. You can change the cycle and syringe entries.

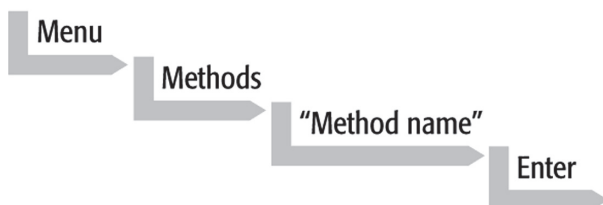
3. If the method is new (that is, added), select and enter a cycle that is appropriate for the application.
4. Select the specific syringe to be used by the method.

**Note** Once you create and save a method, you cannot change the cycle and syringe. To use a different cycle or syringe, you must create a new method.

5. Assign parameter values according to the application requirements.

## Editing and Viewing Methods

You can view and change method parameters (cycle and syringe not included) from the Method menu as follows.

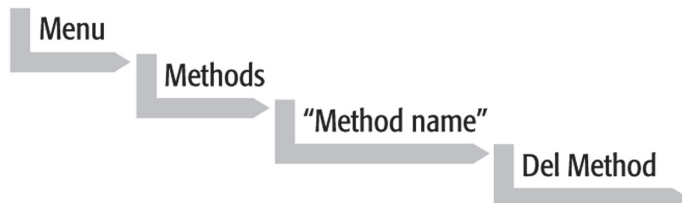


#### ❖ To edit or view a method

1. Scroll to and select the parameter you want to change. Assign the new value and press the ENTER key.
2. Exit the parameter list by pressing either Home function key (F4) to return to the top-level Job Queue menu or the ESC key to return to the previous menu.
3. To view method contents from the job queue displays, select the desired job, press ENTER, and then press the View Method function key.

## Deleting Methods

You can delete methods from the Methods menu, but you cannot delete methods in use by an active job. Complete the following menu selections to delete a method.



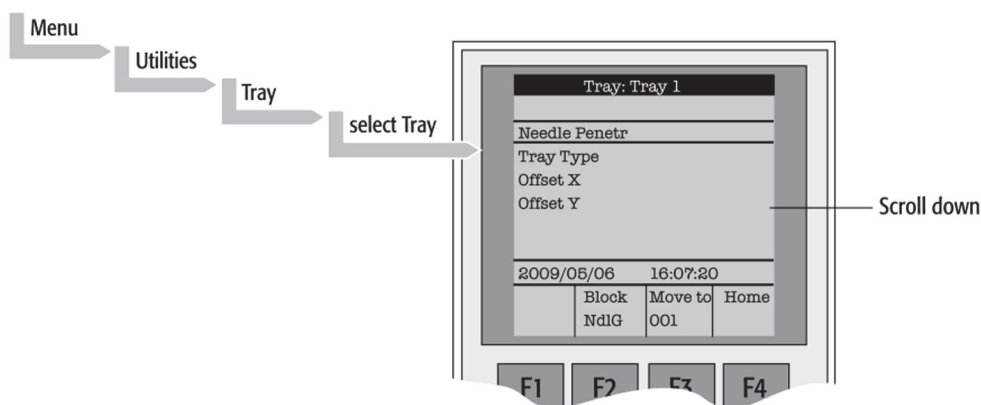
## Job and Job Queue

A job bundles the specified tray with the designated vials (samples) and with the method to run those samples. Another term often used for job in the chromatographic field is *sequence*.

If you prepare more than one job, use the term, *job queue*.

Before activating a job, you must verify that the tray type matches the specified tray and vial size (type) (Figure 51). Perform this step in Utilities.

**Figure 51.** Verify tray with corresponding tray type



Select the corresponding tray type. By activating the function key F3, you can move the injection unit to the first position and to another two corner positions to verify the correct selection of the tray type and teaching position.

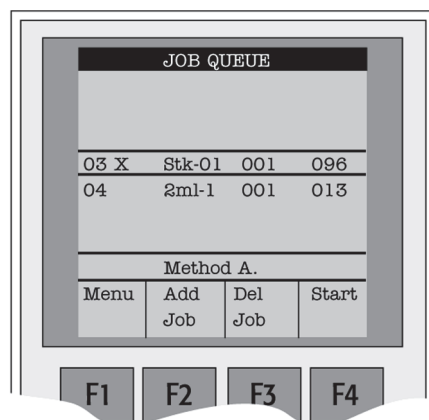
## Building and Starting a Job Queue

### ❖ To build and start a job queue

1. Power on the Accela Open Autosampler.

The JOB QUEUE display opens (Figure 52).

**Figure 52.** JOB QUEUE



2. Load a sample tray onto an available location in a tray holder or stack. Note the corresponding tray name.
3. Add a new job for the tray. Press the Add Job key to bring up the default job.
4. For a tray, select the tray name (for example, stk1-01) that corresponds to the location of the tray you load.
5. Enter the first and last sample number for this job.
6. Select and enter the sample processing method for this job.
7. Press the Home function key (F4) to return to the JOB QUEUE display.
8. To add additional samples that you want to process, repeat steps 2 through 7.
9. If necessary, replace or clean the syringe (liquid versions only), or do both.

Press Menu to see the available options for changing (F1: Change Syringe) and cleaning (F2: Clean Syringe). To completely remove air bubbles, prime the syringe manually.

10. If processing only one job, select the job with the scroll knob. Press the Start key. In the Select Job(s) to Process dialog box, select one of the following options:
  - **All** (entire job queue starting from the top)
  - **Selected** (job selected with the cursor bar)
  - **Resume** (continue with the next job after the one that was canceled)

## Canceling a Job Queue

### ❖ To cancel a job queue

1. In the JOB QUEUE display, press STOP.
2. Select one of the available options:

Continue, Sample, Job, or Job Queue.

- To resume processing with the current sample, select **Continue**.
- If there is a problem with only the current sample, select **Sample**. Processing resumes with the next sample.
- To cancel processing all samples in the current job, select **Job**. Processing resumes with the next job. The canceled job is marked with an X.
- To cancel processing all jobs, select **Job Queue**. The JOB QUEUE display appears. The canceled job is marked with an X.

## Restarting a Canceled Job Queue

### ❖ To restart a canceled job queue

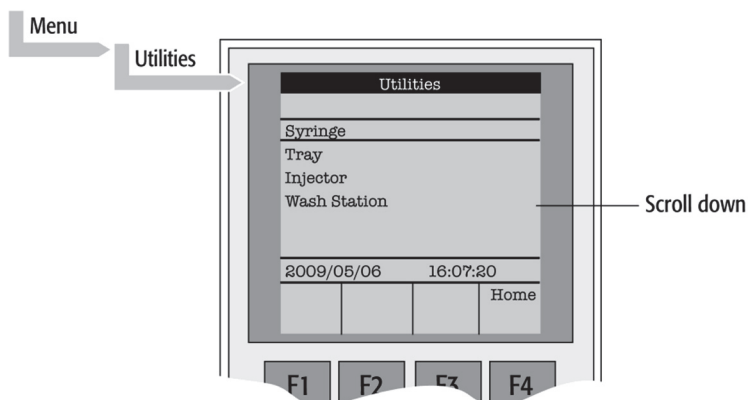
1. In the JOB QUEUE display, press START.
2. Select **Resume**. The job after the last one marked as canceled starts.

## Utility Functions

Utility functions, selectable from the Menu display (Figure 53), provide quick access to checking operations and parameters that you might need to change. These functions are available for the actual syringe, trays, injectors, and wash station. They provide access to key functions without having to set up and execute a method and job.

**Note** If the sample processing cycle uses an item, the method value overwrites the appropriate utility value.

**Figure 53.** Selecting Utilities functions



## Syringe

Table 8 lists the functions available for Syringe by pressing the appropriate function key.

**Table 8.** Utility function keys for Syringe

Function key	Description
F1 Chang Syr	The syringe is moved to a position where the syringe assembly can be completely lowered to ease removal of the syringe adapter. You can then remove the syringe from the adapter and replace it. A prompt is displayed to specify the new syringe. Install the syringe before pressing Enter.
F2 Clean Syr	Use to clean or prime the syringe prior to use. After pressing F2, select either Wash1 or Wash2.
F3 Set Pos	Use to define the Chang Syr position.
F4 Home	The injection unit moves to its Home position and the Job Queue menu displays.

Table 9 lists the items you can select to change the syringe configuration.

**Table 9.** Syringe items (Sheet 1 of 2)

Item	Description
Actual ID	Indicates the identification number (ID) of the currently inserted syringe. If the syringe detection system is set to manual, this message displays: Syringe: No syringe.
Fill Volume	Controls the filling of the syringe. Air bubbles might remain below the plunger after the first pull up. If the plunger moves up and down several times, these air bubbles work out. With this operation you can completely fill the syringe even when using very small sample volumes.



**Table 9.** Syringe items (Sheet 2 of 2)

Item	Description
Fill Strokes	Number of fill strokes. All fill strokes, except the last one, use the selected fill volume. If the selected sample volume is greater than the fill volume, the sample volume is used for all fill strokes. If you select zero, the plunger pulls up only once using the sample volume value.
Pullup Del	Allows you to select a delay time between sample pullup and ejection while filling the syringe. When the plunger reaches the zero position during the fill strokes, the system waits half the Pullup Del time. This allows for an air bubble to float away from the needle tip. This feature is especially useful for removing any air bubble in the syringe and handling viscous fluids.
Fill Speed	Speed of plunger movement used in all syringe filling operations.
Eject Speed	Speed of plunger movement used in all syringe eject operations except sample injection.
Inject Speed	Speed of plunger movement for sample injection. Typically used for Fill Strokes.
Plunger Chnge Pos	Plunger position during the Change Syringe operation. The syringe plunger moves to a position where you can remove and replace the syringe. You may change the value for different types of syringes.

## Tray

Table 10 lists the functions available for Tray by pressing the appropriate function key.

**Table 10.** Utility function keys for Tray

Function key	Description
F2 Block NdlG	Activates needle guide blocking. The option after activation is Rel NdlG, which releases needle guide blocking. Use this function key to test the functionality of the solenoid that blocks the needle guide.
F3 Movto nnn	Serves as a quick check to determine if the <i>x</i> , <i>y</i> , <i>z</i> coordinates are defined correctly for the selected tray. To use this utility, the selected tray, including the sample vials, must be present. After pressing Movto 001, the injection unit moves to sample position No.1. You can repeat this procedure for the last sample position in the first row and the last sample position.
F4 Home	The injection unit moves to its Home position and the Job Queue menu opens.

Table 11 lists the items you can select to change the tray configuration.

**Table 11.** Tray items

Item	Description
Needle Penetr	Needle penetration depth into the sample vial. You can change the needle penetration depth for the selected tray by entering the value that you want.
Tray Type	Shows the tray type. If the tray enables the use of different tray types, you can make the change at this position.
Tray Offset X	Use to correct the ideal $x$ position of Position 1 if necessary.
Tray Offset Y	Use to correct the ideal $y$ position of Position 1 if necessary.
Tray Offset Z	Use to correct the ideal $z$ position of Position 1 if necessary.
dxRow	Corrects any inclination of a tray (plate) in the $x$ axis of a row.
dyRow	Corrects any inclination of a tray (plate) in the $y$ axis of a row.
dzRow	Corrects any inclination of a tray (plate) in the $z$ axis of a row.
dxCol	Corrects any inclination of a tray (plate) in the $x$ axis of a column.
dyCol	Corrects any inclination of a tray (plate) in the $y$ axis of a column.
dzCol	Corrects any inclination of a tray (plate) in the $z$ axis of a column.

## Injector

Table 12 lists the functions available for Injector by pressing the appropriate function key.

**Table 12.** Utility function keys for Injector

Function key	Description
F3 Movto Inj	The injection unit moves to the selected injector position. With this function, for example, the injectors GC-Inj1, Waste, Waste2, and Flush can be accessed. By selecting the Needle Penetr parameter on the same screen, you can check or change the injector needle penetration value.
F4 HOME	The injection unit moves to its HOME position and the Job Queue menu is displayed.

Table 13 lists the item you can select to change the injector configuration.

**Table 13.** Utility item

Item	Description
Needle Penetr	By selecting the Needle Penetr parameter, you can check or change the injector needle penetration value. To ensure reproducible sample injections and minimize carryover, make sure that the needle penetration depth is accurately set.

## Wash Station

Table 14 lists the functions available for Wash Station by pressing the appropriate function key.

**Table 14.** Utility function keys for Wash Station

Function key	Description
F3 Movto Wash	The injection unit moves to the selected wash station port. By selecting the Needle Penetr parameter on the same screen, you can check or change the wash station needle penetration value.
F4 HOME	The injection unit moves to its HOME position and the Job Queue menu is displayed.

Table 15 lists the items you can select to change the wash station configuration.

**Table 15.** Wash station items

Item	Description
Needle Penetr	By selecting the Needle Penetr parameter, you can check or change the wash station needle penetration value.
Rinse Time	If rinse time is activated (value > 0), the solenoid of a fast or active wash station opens for the specified time after the syringe needle has been removed from the wash port (after completion of the syringe wash cycle).  The solvent flows into the wash port without the restriction of the needle; be aware of higher solvent consumption.

## Vial

Table 16 lists the functions available for a vial by pressing the appropriate function key.

**Table 16.** Utility function keys for a vial

Function key	Description
F3 Movto Vial	The injection unit moves to the selected vial type. By selecting the Needle Penetr parameter, you can check or change the wash station needle penetration value.
F4 HOME	The injection unit moves to its HOME position and the Job Queue menu is displayed.

Available Vial types:

- Standard: Vial in standard wash station for gas chromatography (front position).
- Fiber Exp: Position to block needle guide. Mainly used for SPME application. Refer to the SPME User Manual.

Table 17 lists the item you can select to change the vial configuration.

**Table 17.** Vial item

Item	Description
Needle Penetr	By selecting the Needle Penetr parameter, you can check or change the wash station needle penetration value.

## Dilutors

Table 18 lists the functions available for dilutors by pressing the appropriate function key.

**Table 18.** Utility function keys for dilutors

Function key	Description
F1 Prime	Primes the dilutor syringe with solvent.
F2 Chang DSyr	Moves the plunger of the dilutor syringe to standby position to allow easy access to change the syringe.
F4 HOME	The injection unit moves to its HOME position and the Job Queue menu is displayed.

Table 19 lists the items you can select to change the dilutor configuration.

**Table 19.** Dilutor items

Item	Description
Syringe	Indicates the dilutor side-port syringe inserted in the z axis. Use this function to select another syringe size.
Syr Dilut Pos	Moves the plunger of the side port syringe up by the specified distance. This allows adjustment of the plunger tip of the side port syringe exactly above the lower side port (solvent inlet). You must fine-tune after changing a side port syringe to allow unrestricted solvent flow.
Dilutor Syr	Indicates dilutor syringe is installed. If you change the syringe size, this item must be adapted accordingly. This function coordinates the syringe-specific dimensions.
Prime Volume	Specifies the volume to prime the dilutor syringe can be specified. The allowed range can be fourfold higher than the actual syringe volume. This allows filling and emptying a dilutor syringe more often than just once.
Pullup Delay	Permits selecting a delay time between solvent filling and ejection while filling (or priming) the syringe.
Fill Speed	Speed of plunger movement of dilutor syringe used in all syringe filling operations.
Eject Speed	Speed of dilutor syringe plunger movement used in all dilutor syringe eject operations.
Eject Delay	Permits selecting a delay time between solvent ejection and filling while filling (or priming) the syringe.

## Tools

**Note** The Object class Tools is available with PAL Firmware level 3.0.x, 4.1.x, or later.

Table 20 lists the functions available for tools by pressing the appropriate function key.

**Table 20.** Utility function keys for tools

Function key	Description
F1 Chk Offs	Use Check Offset to verify the offset from the syringe needle tip to the tool tip. This function is mainly used for the MALDI tool; it is not active for the MHE tool.
F4 HOME	The injection unit moves to its HOME position and the Job Queue menu is displayed.

Table 21 lists the items you can select to change the tools configuration.

**Table 21.** Tools items

Item	Description
Teach Point	The reference point at which to teach the Object (Tool; MHETool) is selectable. The MHETool needs no extra position. The Parking Station is the fix point. Select None.
ToolOffset X	Corrects the ideal <i>x</i> position of the MHETool if necessary.
ToolOffset Y	Corrects the ideal <i>y</i> position of the MHETool if necessary.
ToolOffset Z	Corrects the ideal <i>z</i> position of the MHETool if necessary. You must have a ToolOffsetZ of -48.0 mm to position the MHETool on the sample vial.

## Logfile

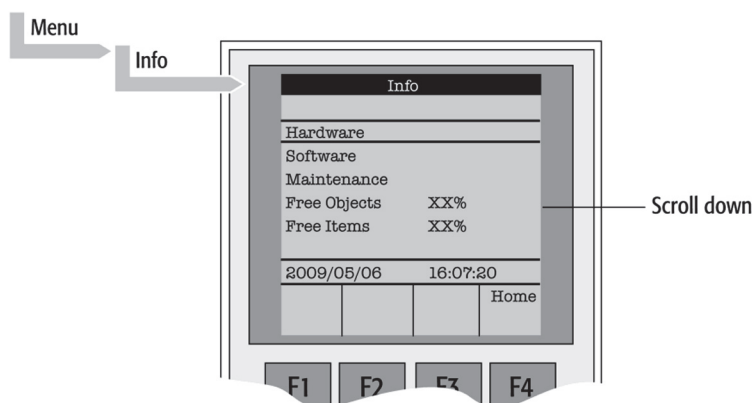
You cannot directly read the PAL logfile on the terminal display. However, the F3 function key does support printouts. You must have a serial printer connected to port SER2 for the PAL System and to port SER1 for PAL-xt system. If a serial printer is not available, use a serial or parallel converter.

## Info Functions

Info functions, selectable from the Menu screen (Figure 54), provide quick access to viewable information. These Info functions are available for the hardware, software, maintenance, free objects, and free items.

All items with the exception of maintenance are read-only.

**Figure 54.** Selecting Info functions



## Hardware

Table 22 lists Info functions items for hardware.

**Table 22.** Info functions items for hardware

Item	Description
CPU SNo	The serial number (S/N) of the PCB APR CPU is displayed.
CPU ID	Version number of the PCB APR CPU.
MOTIO ID	Version number of the PCB APR CPU.

## Software

Table 23 lists Info functions items for software.

**Table 23.** Info functions items for software

Item	Description
PAL Firmware	Firmware level of PAL System
Head Firmware	Version level of the firmware PAL Injection Unit
Terminal FW	Version level of the firmware PAL Terminal
Altera Firmware	Version level of the Altera component firmware

## Maintenance

Table 24 lists Info functions items for maintenance.

**Table 24.** Info functions items for maintenance (Sheet 1 of 2)

Item	Description
PlgStrokeCnt	<p>A counter for syringe plunger movements. The actual number of strokes is displayed. If the counter reaches the set limit (PlgStrikeLim), a warning is displayed at the next start of a job (run). The system continues but signals the user to verify syringe conditions. You can restart the counter by setting it back to zero.</p> <p>A Syringe has only one counter. If you change syringe types, the system continues to count as if it were the same type.</p>
PlgStrokeLim	Sets an upper limit for the syringe plunger strokes.

**Table 24.** Info functions items for maintenance (Sheet 2 of 2)

Item	Description
Inject Count	Monitors the number of injections. The number of injection valve switches is a helpful tool for deciding on replacement parts for the injector system. The actual number of injections (valve switches) is displayed. If the counter reaches the set limit (Inject Limit), a warning appears at the next start of a job (run) (Inject Limit). The same counter is used for injector penetrations with a GC technique.
Inject Limit	Sets an upper limit for the number of injections.

**Note** Counters for the plunger movement and injector penetrations are available with PAL Firmware level 2.5.x or later.

## Free Objects/Free Items

In addition to the core software, the PAL Firmware contains data for firmware objects. There are different classes of Objects, such as Syringes, Trays, Tray Holders, and so on. Each class of objects contains items. The items contain the actual data, such as  $x$ ,  $y$ ,  $z$  positions.

The data are stored in a flash memory backed up by a battery.

To optimize RAM and Flash memory use, the software has a reserved section of memory for each of the Objects and Object Items.

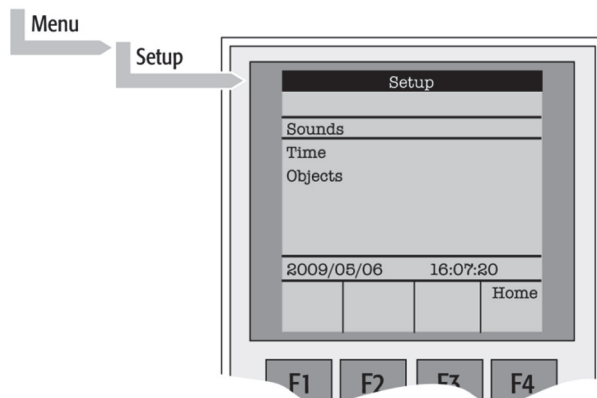
The percentage shown in the Info section indicates how much of the reserved software space is still available.



## Setup Functions

The Setup functions, selectable from the Menu screen (Figure 55), allow access to various functions for the PAL System. The Sound, Time, and Objects are basic functions that you use at installation or if changes have been made over time.

**Figure 55.** Selecting Setup functions



## Sounds

Table 25 lists Setup functions for Sounds.

**Table 25.** Setup functions for Sounds

Item	Description
Message Box	A specific dual beep tone signals that a message box for user intervention appears on the screen. You can turn this beep signal on or off.
Warn Move	Sounds a beep at the start of the PAL movement. Keep this function turned on for safety reasons.
End Cycle	Sounds a beep at the end of a cycle. Select as desired.
End Job	Sounds a beep at the end of a job. Select as desired.

## Time

Table 26 lists Setup functions for Time.

**Table 26.** Setup functions for Time (Sheet 1 of 2)

Item	Description
Year	Sets the year for the PAL internal clock.
Month	Sets the month for the PAL internal clock.
Day of Month	Sets the day for the PAL internal clock.

**Table 26.** Setup functions for Time (Sheet 2 of 2)

Item	Description
Hours	Sets the hours for the PAL internal clock.
Minutes	Sets the minutes for the PAL internal clock.
Seconds	Sets the seconds for the PAL internal clock.

After setting or resetting the date and time, use the F1 function key (F1 Set Time) to save your changes.

## Objects

You can directly access common functions for the various PAL Firmware Object classes. For a detailed listing, see “[PAL Firmware Overview](#)” on [page 136](#).

## Communicating with the Computer

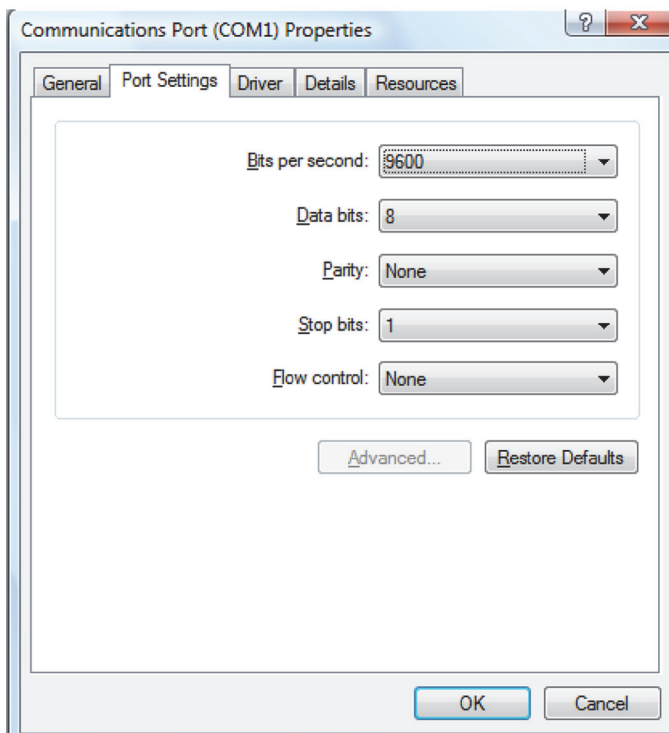
This section describes how the Accela Open Autosampler communicates with a computer.

The PAL System provides a serial communication protocol with a computer. The Accela Open Autosampler system requires a dedicated APR Control-xt board and PAL Firmware level 4.1.x or later.

The COM port settings are normally defined in the Windows Device Manager (choose Control Panel > Hardware and Sound [or Device Manager if set to the icons view] > Ports).

[Figure 56](#) shows the standard settings for the serial communication. Be aware that the application (for example, PAL Loader software) does actively set the communication parameters, and the baud rate is set for optimized use.

**Figure 56.** COM Port Settings



A quick way to check whether serial communication with the PAL System can be established is to use the PAL Loader software. For further information, see “PAL Loader” on [page 155](#).



# Operating the Temperature Controllable Stack Coolers

This chapter describes how to set the temperature, temperature control, and alarm. It also discusses temperature stability and condensation build-up.

## Contents

- [Temperature Setting, Control, and Alarm](#)
- [Temperature Stability](#)
- [Condensation Build-Up](#)

## Temperature Setting, Control, and Alarm

This section contains information about the temperature setting, temperature control, and temperature alarm.

### Temperature Setting

Turning the power on (I) at the power supply module puts the control unit into self-check mode. The set value for the temperature appears after a few seconds.

The stack/tray cooler temperature is preset by the factory to 10 °C.

#### ❖ To change the value

1. Press **P** briefly (less than two seconds) until SP1 appears.
2. Use the increment/decrement keys to select the desired value.
3. Press **P** again.

## 4 Operating the Temperature Controllable Stack Coolers

Temperature Setting, Control, and Alarm

**Note** Pressing the P button for two seconds or longer will access the programming level. You can access values, such as OPEr, ConF, OFF, rEG, tune and OPLO, in the menu by selecting the increment/decrement keys. The user is blocked from all of these menu items. Leaving the unit untouched for approximately two minutes returns the display to the current temperature reading information. Turning the device off/on reinstates the temperature display.

Factory programming blocks the U key.

There are indicators below the display digits that show which relay is in use, OUT1 or OUT2. The arrows indicate if the unit is heating or cooling.

◀ heating      — temperature within range of set point      ▶ cooling

## Temperature Control

As described in the Specifications section, the cooling/heating capacity of a temperature control device based on the Peltier technique is always related to room temperature. Only a limited maximum temperature difference from ambient temperature can be achieved. PAL stack/tray cooler modules should reach a minimum difference of 14 °C ( $\Delta T$ ).

The controller provides the following temperature setting range:

- Stack cooler: 1 to 45 °C
- Tray cooler: 1 to 70 °C

This wider control range, which is greater than the specification limits (4 to 40 °C and 4 to 70 °C, respectively), provides cooling or heating when the ambient temperature is too high or too low to easily reach the desired temperature. Heating to a higher temperature could be important for metabolism or kinetics studies at the human body temperature level of 37 °C.

The cooling option is usually necessary to protect the sample from ambient temperature during the analytical run time. Stack/tray cooler devices are rarely used to store a sample at a given temperature for a longer time period.

The measured and displayed temperature represents the temperature in the compartment and not in the sample liquid. To check the temperature control, an independent temperature probe is inserted from the front by removing one plastic screw from the front cover. Position the probe in the middle of the drawer where the tray separator projects up. Tape the probe to the metal plate of the drawer.

To reach 4 °C in the analytical solution, program a set value lower than 4 °C. The material used for the sample may be glass, polypropylene, polyethylene, or similar polymer products. All of these have excellent insulating properties.

If the sample has to be cooled as low as 4 °C, cool the sample tray first in a refrigerator before the tray is placed into a stack/tray cooler module. This shortens cool-down time considerably.

## Temperature Alarm

An internal alarm point of 55 °C protects the upper limit of a stack cooler module; it is 72 °C for a tray cooler. Reaching this point in a single fault situation causes the temperature control unit to automatically force cooling to the fullest extent possible. (The alarm is not audible.) At the same time a corresponding Error message is displayed. For explanations of these messages, see “Cooled Stack Power Supply Error Messages” on page 90.

A fuse to prevent overheating is built in for unattended automated runs.

The Peltier element turns off automatically at  $72 \pm 5$  °C.

Only an authorized representative of Thermo Fisher Scientific can replace a damaged over-temperature fuse.

In case of error, do the following systematically:

1. Check the corresponding error message (“Cooled Stack Power Supply Error Messages” on page 90).
2. Verify whether the appropriate power supply is being used for the stack cooler and tray cooler.
3. Turn off the power supply, wait a few minutes, and turn it on again.

Ensure that the alarm is off. If the alarm is still on, observe the corresponding error message.

## Temperature Stability

The stack/tray cooler is basically a thermostatted sample tray that keeps the analytical sample below or above ambient temperature.

Thermo Fisher Scientific recommends the following guidelines:

- Switch on the stack/tray cooler at least 30 minutes before the analytical routine run is to be started at 10 °C or 75 minutes, if operated at 4 °C.
- For high-throughput analysis—typical cycle time of 60 seconds or less per analysis—leave the drawers open between injections.
- For longer cycle times, use the Accela Open Autosampler option to close the drawer after each sampling.

You can change the Restore mode to keep the drawer open or closed during a run.

**Note** The preset mode for the stack cooler in combination with the PAL System closes the drawers after sampling during an analytical run.

❖ **To change the Restore mode**

Complete the following steps starting at the top-level command from the control terminal:

- a. Choose **Menu > F3-Setup > Objects > Trayholders > Cstack > Restore Mode**.
- b. Change the default value from Sample to **Auto**.
- c. Press HOME.

## Condensation Build-Up

This section contains information about condensation build-up in stack coolers and tray coolers.

### Condensation Build-Up in Accela Open Autosampler Stack Coolers

Condensation build-up is directly related to the temperature and relative humidity/temperature in ambient air (dew point). Long-term tests showed very little build-up of condensation in an environment of relative humidity up to 60 percent and  $22 \pm 2$  °C ambient temperature.

Condensation at the back of the compartments is channeled to the drain outlet labeled Condensation Drain at the lower back portion of the stack cooler. The outlet is plugged with a paper filter. This helps to evaporate the collected water in the drain line using the excess heat from the Peltier element. A drain line from the outlet to a reservoir bottle is not necessary under normal conditions.



**CAUTION** If a flush gas is used as described below, you are responsible for ensuring that a two-stage safety pressure regulator device is installed between the gas supply and the stack cooler.

Do not use any flammable or explosive gas such as hydrogen.

If you operate the stack cooler under severe conditions, you can use a flow of dry and clean (oil-free) air or nitrogen to dry the compartments continually. Connect the corresponding gas line to the Swagelok fitting (1/8 in.) at the back of the stack cooler DW labeled Flush Gas. A flow of approximately 300 to 400 mL/min is required to keep the compartments moisture-free.

If acidic vapor phase is anticipated due to the application, use the same gas line connection to flush a stream of inert gas, such as nitrogen or helium, into the compartments. A stream of a few mL/min can help to prevent corrosion of the rolls.



Thermo Fisher Scientific recommends checking regularly for condensation build-up. It is good practice to clean the inside of the stack cooler when changing the analytical samples, dry out the stack cooler at ambient temperature at regular intervals, and open the drawers 1–5 cm (0.5–2.0 in.) for air circulation. The compartment drying frequency depends greatly on surrounding conditions.

## Condensation Build-Up in PAL Tray Coolers

Condensation build-up is directly related to the temperature and relative humidity/temperature in ambient air (dew point). Long term tests showed little build-up of condensation in an environment of relative humidity up to 60 percent and  $22 \pm 2$  °C ambient temperature.

A built-in drain line, as installed in stack coolers, is not provided in a tray cooler.

It is also good practice to install the Plexiglass lid whenever possible to avoid condensation build-up. You might not be able to install a lid for certain applications, for example, if a vial has to be transported to an Agitator or a Barcode Reader. In certain situations, the height of the vial is the limiting factor for installing a lid. Operating a tray cooler under less than optimal conditions requires some on-site improvisation. A simple foil or lid with cut-outs for the vials will limit condensation build-up dramatically.

Thermo Fisher Scientific recommends checking regularly for condensation build-up. Clean the inside of the tray cooler when changing the analytical samples, and dry out the tray cooler at ambient temperature at regular intervals.

**Note** For certain applications, a change from a tray cooler to Stack 2DW Cooler might be indicated. As long as deep well, microtiter plates, or 2 mL vials (TrayType VT54) are being handled, this change is positive with respect to condensation build-up. It would also double the capacity of the trays.



## Maintenance

This chapter describes the maintenance procedures that can help you operate the Accela Open Autosampler under optimal conditions.

To ensure accuracy and precision of the Accela Open Autosampler. Follow the maintenance procedures at the suggested intervals in [Table 27](#).

If you use the system extensively (for example, nights and weekends), or if you use corrosive solvents, you might need to perform the maintenance procedure more frequently. See [Table 28](#).

**Table 27.** Routine maintenance for Accela Open Autosampler (Sheet 1 of 2)

Maintenance step	Interval
Clean the outside of the instrument. Use only a soft, lint-free cloth dampened with mild soap and water.	Weekly or as needed
Clean instrument, syringe, and surfaces.	Weekly or as needed
Replace the valve needle seal.	Check the tightness of the needle in the needle seal.
Replace the valve rotor.	Annually or more often, depending on the throughput and quality of sample solution and mobile phase
Clean the outside of the instrument. Prime the entire DLW system by using the dedicated macro (cycle).	Daily or as needed (required before routine running)
Replace the DLW syringe plunger.	The syringe plunger for a gas tight syringe (polymer tip) has to be replaced on a regular basis. The interval is highly dependent on the application, throughput, and quality of the sample solution (particles, and so on) washing solvent. Check the tightness of the plunger on a regular basis to gain experience with the application.

**Table 27.** Routine maintenance for Accela Open Autosampler (Sheet 2 of 2)

Maintenance step	Interval
Replace the DLW syringe needle. Observe the correct placement of the Teflon seal.	Only as needed (for example, if bent)
Check all connections for tightness.	Weekly or as needed
Replace the DLW holding loop.	Recommended once a year as preventative maintenance
Check the spring-loaded needle holder assembly for functionality. Check the spring compression in particular.	Weekly or as needed
Wash and waste solvent reservoirs: check flow restriction of frits.	Weekly or as needed
Quality of solvents, biological growth	
Check the functionality of the DLW actuator/solenoid.	Weekly or as needed
	Change once a year as preventative maintenance.

Specific kits for the GC or the HPLC techniques are available. The kits include parts that you must change annually.

## Cooled Stack Maintenance

If you use the system extensively (for example, nights and weekends), or if you use aggressive solvents, you might need to perform maintenance procedures more frequently.

**Table 28.** Stack/tray cooler maintenance (Sheet 1 of 2)

Maintenance step	Interval
Clean the outside of the instrument. Use only a soft, lint-free cloth dampened with mild soap and water.	Weekly or as needed
Clean and dry the inside of the instrument if necessary, including compartments, rolls, and rails. Use only a soft, lint-free cloth.	Daily or as needed

**Table 28.** Stack/tray cooler maintenance (Sheet 2 of 2)

Maintenance step	Interval
Clean and grease the rolls and the rails <sup>a</sup> .	<p>Once a year without acidic vapor exposure</p> <p>If acids are used for the sample solution and the stack cooler is exposed to the vapor, clean more frequently, as required.</p> <p>A positive stream of inert gas, such as nitrogen, dried and clean air, or helium, can prevent corrosion. See “<a href="#">Condensation Build-Up</a>” on page 78.</p>
Verify the position and condition of the stopper and holding magnet of the drawer or drawers.	Monthly or as needed

<sup>a</sup> Only valid for stack cooler modules

**Note** Recommended Grease and Supplier:

- Grease (PN: CH-952961): approx. 8 g grease and brush
- Grease Temperature Range: –130 to +150/+200 °C
- Supplier: Thermo Fisher Scientific

**Note** There are no operator-serviceable or replaceable parts inside the power supplies or the PAL System. In case of failure, contact a Thermo Fisher Scientific representative.

**Note** In the Power Supply (PN CH-952995), a fan is installed in the housing to cool the electronic components. Fan cooling prolongs the life of the electronic components.



**CAUTION** Check the fan regularly. If the fan fails, the power supply still operates and can pass safety tests, even when the temperature is dangerously high and might damage the system electronics.



## Troubleshooting

This chapter describes the troubleshooting procedures based on different error messages. It also includes some special considerations for troubleshooting and standard chromatographic tests.

### Contents

- [Troubleshooting Using Observed Symptoms or Error Messages](#)
- [Cooled Stack Power Supply Error Messages](#)
- [Troubleshooting Considerations](#)
- [Standard Chromatographic Tests](#)

## Troubleshooting Using Observed Symptoms or Error Messages

Table 29 lists the symptoms or error messages and how to troubleshoot them.

**Table 29.** Troubleshooting using symptoms and error messages (Sheet 1 of 5)

Symptom or error message	Possible cause	Recommended action
Detector signal not detected or very low	Clogged syringe	Remove syringe and aspirate/dispense liquid manually. Clean syringe.
	Bent needle	Inspect and change syringe if necessary.
	No sample liquid injected	Check and adjust Needle Penetration into sample vial if necessary.
	Sample volume too low	Increase sample volume.
	Improperly installed valve needle guide and/or needle seal	Check valve needle guide and seal.
	Valve ports not plumbed correctly to the pump, the detection system, or both	Check plumbing connections.
	Wrong valve type specified	Check valve type by choosing <b>Menu &gt; Setup &gt; Objects &gt; Injectors &gt; LCV1v1 &gt; Valve.</b>

## 6 Troubleshooting

Troubleshooting Using Observed Symptoms or Error Messages

**Table 29.** Troubleshooting using symptoms and error messages (Sheet 2 of 5)

Symptom or error message	Possible cause	Recommended action
Flow of mobile phase and/or sample from the wrong ports on LC injection valve	The valve rotor is reversed (180 degrees out of alignment).	Remove the rotor and reinstall in the correct position. For a Cheminert valve, check for marking points. For a W-type valve, look for a letter (for example, H) stamped on the rotor. The letter must be on the lower half of the rotor.
Detector signal not detected or very low	Clogged syringe, needle, or holding loop	Remove syringe and aspirate/dispense liquid manually. Clean syringe, needle/holding loop, or both of these.
	Bent needle	Inspect and change syringe if necessary.
	No sample liquid injected	Check and adjust Needle Penetration into sample vial if necessary.
	Sample volume too low	Increase sample volume.
	Improperly installed valve needle guide, needle seal, or both	Check valve needle guide and seal.
Sample backing up on the valve needle guide	Valve ports not plumbed correctly to the pump, detection system, or both.	Check plumbing connections.
	Wrong valve type specified	Check valve type by choosing <b>Menu &gt; Setup &gt; Objects &gt; Injectors &gt; LCVlv1 &gt; Valve.</b>
	Syringe OD too small	Check the syringe needle for the correct gauge. Gauge 22 = 0.72 mm OD
	Leaking valve needle seal	Change valve needle seal.
	Needle penetration depth set incorrectly for the injection valve	Adjust the injection valve needle penetration.
	The syringe plunger speed set too high, resulting in excessive pressure in inlet	Reduce inject speed in method.



**Table 29.** Troubleshooting using symptoms and error messages (Sheet 3 of 5)

Symptom or error message	Possible cause	Recommended action
Syringe not filling properly	DLW Actuator/Solenoid not functioning	Load PAL Firmware level smaller than 4.0 or PAL Firmware Object for DLW option onto the system.
		Solvent path is blocked. Prime the system or inspect it to identify the reason for clogging.
		Air bubbles are trapped in the system, caused by loose connections or tubing not cut square.
	Worn out syringe plunger tip	Replace syringe plunger. Check if glass barrel is scratched (damaged). If in doubt, replace the entire syringe at once.
	DLW Actuator Control PCB defect	If possible, verify the DLW Actuator Holder assembly on another PAL System. If defect is definite, replace PCB.
	DLW Flow Diverter	Check the Flow Diverter: <ul style="list-style-type: none"> <li>• Is the Flow Diverter inserted? Inserted in full up position?</li> <li>• Is the solvent flowing freely in the tube?</li> </ul>
	Fitting replaceable needle	Check the fixation and seal of the replaceable syringe needle: <ul style="list-style-type: none"> <li>• Is the retaining nut tight?</li> <li>• Is the PTFE seal installed?</li> <li>• Is the needle entry free of debris, not clogged?</li> <li>• Is the correct needle diameter (gauge 22) used?</li> </ul>

## 6 Troubleshooting

Troubleshooting Using Observed Symptoms or Error Messages

**Table 29.** Troubleshooting using symptoms and error messages (Sheet 4 of 5)

Symptom or error message	Possible cause	Recommended action
Sample peaks/responses not reproducible	Dirty syringe or Holding Loop	Increase Pst Cln Slv1/Slv2 values in method. Use PreClnSlv1/Slv2 and PreClnSpl.
	Syringe pressure differences	Increase Pullup Delay value.
	Vacuum created in sample vial	Reduce sample volume in sample vial. Use setting under <b>F3-Setup &gt; System &gt; PrePressureVial</b> .
	Method parameters	Check recommended Method parameters in “PAL Firmware Overview” on page 136. First verify these critical parameters: <ul style="list-style-type: none"><li>• Fill speed</li><li>• Pullup delay</li><li>• Injection Speed</li><li>• Post Inj Delay</li></ul>
Excessive carryover between samples	Loose, unstable, or wrong connection	Check all connections within the DLW system, such as needle, Holding Loop, injection loop connections, and so on.
	Inappropriate wash solvent or solvents	Use appropriate wash solvents.
	Waste tubing ID too small at injection valve, causing waste liquid to be pulled back, by capillary action, into rotor groove	Replace the waste tubing with larger ID tubing.
	Air gaps not formed. Liquids of different type blending into each other.	Verify the DLW method parameter for Airgap Volume.  Prime entire system.

**Table 29.** Troubleshooting using symptoms and error messages (Sheet 5 of 5)

Symptom or error message	Possible cause	Recommended action
Excessive carryover between samples	Dirty needle, holding loop, and/or valve injection port	Increase Pst Inj Slv1/Slv2 values in method (cleaning time).  Use PreClnSlv1/Slv2 and PreClnSpl.
	Inappropriate wash solvent or solvents	Use appropriate wash solvents.
	Waste tubing ID too small at injection valve, causing waste liquid to be pulled back, by capillary action, into rotor groove.	Replace the waste tubing with larger ID tubing.
	Damaged or grooved valve rotor	Replace valve rotor.
	Leaking valve needle seal	Change valve needle seal.
	Inappropriate wash solvent composition	Use appropriate wash solvent. Also observe the order of use for biological samples. First wash is always an aqueous solution followed by organic solvents.
	Tube of holding loop at end not cut square or nut/ferrule fixation not correct.	Check the connections from the holding loop: <ul style="list-style-type: none"> <li>• Is the pilot distance appropriate?</li> <li>• Is the tube end damaged?</li> <li>• Is the tube end cut absolutely square?</li> <li>• Are the right nuts and ferrules used?</li> </ul>
	DLW Actuator/Solenoid questionable or defective	Verify the functionality of the DLW Actuator.  Is solvent flowing when active? Use command <b>Utilities &gt; Wash Station</b> to test.
	Blocked solvent frits in wash solvent reservoir	Clean the frits in an ultrasonic bath with an appropriate solvent. If flow is not ensured after cleaning, replace frits.  Replace wash solvent and clean reservoir bottle and tubing carefully.

## Cooled Stack Power Supply Error Messages

The temperature control unit for the stack cooler can display feedback messages in the form of error messages. See [Table 30](#).

**Table 30.** Stack cooler controller error messages

Message	Description	Reason	Check/Solution
-----	Error probe	DC cable disconnected or in short circuit	Check connection between power supply and Stack/TrayCooler.
		Connection mistake	Check signal probe (for example: NTC measures the resistance: 25 °C = 10 kOhm).
		Damaged Probe	
uuuu	Under range	Measured variable below the sensor limit	Check the set point for the desired temperature.
oooo	Over range	Measured variable above the sensor limit	Turn the unit off/on. Initialize the Temp. Control Unit to normal level.
			Check the set point for the desired temperature.
LbA	Loop break alarm	Interruption of loop break alarm	Check the set point for the desired temperature.
			Turn the unit on/off. Initialize the Temp. Control Unit to normal level.
ErEP	EPROM	Error in Eprom	Press <b>P</b> .

Other messages, such as ErAT and noAT, are suppressed and are not of use in this application.

**Note** The Temperature Control Unit for the stack/tray cooler does not allow the user to enter different levels of programming functions. The program is factory set, downloaded to each Temperature Control Unit at the factory. A power supply programmed for the PAL Stack Cooler cannot be used for a PAL tray cooler.

## Troubleshooting Considerations

This section contains information about special troubleshooting considerations, including tubing connections and solvent delivery pumps not priming or not functioning.

### Tubing Connections

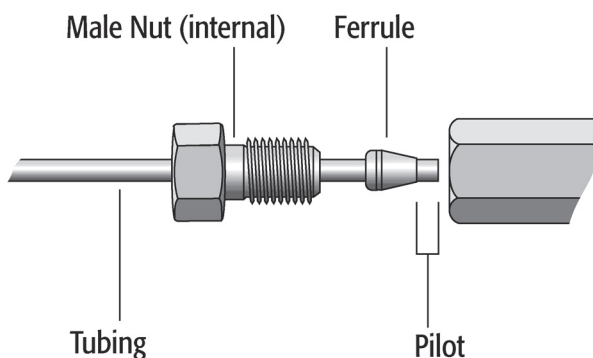
You must pay special attention to all tubing connections, including these common mistakes in routine practice:

- Nut and ferrule not connected leak tight (see [Nut and Ferrule Tightening](#))
- Incorrect pilot distance (see [“Incorrect Pilot Distance”](#) on page 92)

- Tube not cut square (see “[Tube Not Cut Square](#)” on page 93)
- Wrong type of ferrule or nut used (see “[Wrong Nuts or Ferrules Used](#)” on page 94)
- Trapped air bubbles (see “[Avoiding Trapped Air Bubbles](#)” on page 95)

Figure 57 provides an illustration of the terms used for tubing connections.

**Figure 57.** Terms for tubing connections



Such mistakes lead to these issues:

- Dead volume
- Peak deformation or peak splitting
- Carryover effects
- Decreased resolution of the tubing between column and detector

You can prevent issues with tubing connections by observing the following basic rules.

## Nut and Ferrule Tightening

### ❖ To tighten a nut and ferrule stepwise

1. The first important rule when tightening a nut and ferrule is to proceed stepwise.
  - a. Tighten the nut by hand as much as possible.
  - b. Continue to tighten using a wrench until you feel resistance. Add one or two quarter turns to reach the sealing point.
  - c. Open the connection and remove the assembly to inspect the pilot distance. See [Figure 57](#).
  - d. Install the assembly again and, first, manually tighten the nut. Then, use the wrench to reach the same sealing or resistance point as reached in step 2 and add one quarter turn extra for the final seal.
  - e. Check the seal when liquid is pumped through the system. In case a leak is observed, tighten once more by another quarter turn. Always move from a leaking to a tight seal; never move backward from an overtightened seal.

Do not overtighten the nut/ferrule. If a connection proves to be leaky, use another quarter turn to tighten. Step-by-step tightening is the correct approach. Overtightening even once will damage the seat, and the next connection can only be sealed using force. If the seat is damaged, the ferrule will likely stick and require force or special tools (such as a drill) to remove it.

If you use different materials, such as a polymer tubing and nut, together with a stainless steel ferrule, then overtightening is always a danger. The ferrule will bite into the polymer tube and block the flow.

When installing a tube with a narrow bend, typically a loop, fix one connection first; do not simultaneously insert the second nut into the female counterpart. After tightening open the connection as described above, start preparing the second connection without inserting the first nut. Open the second connection again and check the pilot distance. The last step is inserting the complete tubing with the two connections and making a final seal as described earlier.

This stepwise installation of a bended tube is mandatory for PEEKsil tubing (fused silica tubing sheathed with PEEK polymer).

2. The second rule to observe when installing nuts and ferrules is to never reuse them for any other connection.

Do not install a tube for any other application with a nut/ferrule that is set for a specific purpose. Cut a portion of the tube end until straight tubing is reached and remove the ferrule.

This rule is valid for a loop as well. Never reuse a loop for another valve or reuse a valve after exchanging the stator.

If a fingertight nut has been installed, replace it as well. If a stainless steel nut has been used, ensure that you use the correct type.

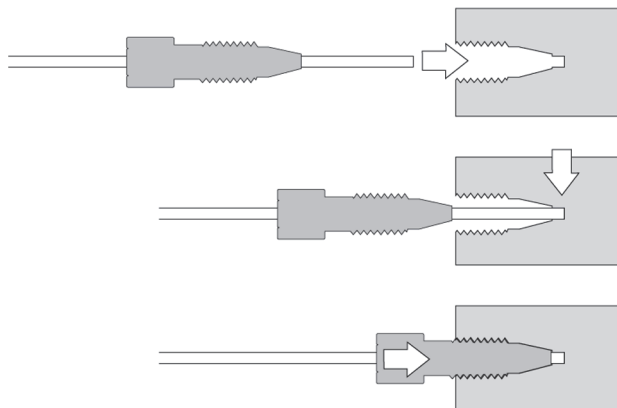
Newly installed fingertight ferrules can hold a backpressure of up to 200 bar.

## **Incorrect Pilot Distance**

When making a new tube connection, be sure to follow these guidelines:

1. Slip the nut and ferrule over the tube. Make sure the end of the tube sticks out approximately 50 mm (2 in.).
2. Press the tube end firmly into the female counterpart for the connection.
3. Move the nut and ferrule together into the female counterpart for the connection (Figure 58). Make sure to maintain pressure on the tube to avoid the tube slipping backwards, out of position.
4. Tighten the nut as described in the previous section.

**Figure 58.** Pressing the tube into place



Failure to observe this procedure will lead to an incorrect pilot distance, which will result in actual dead volume, not delay volume.

## Tube Not Cut Square

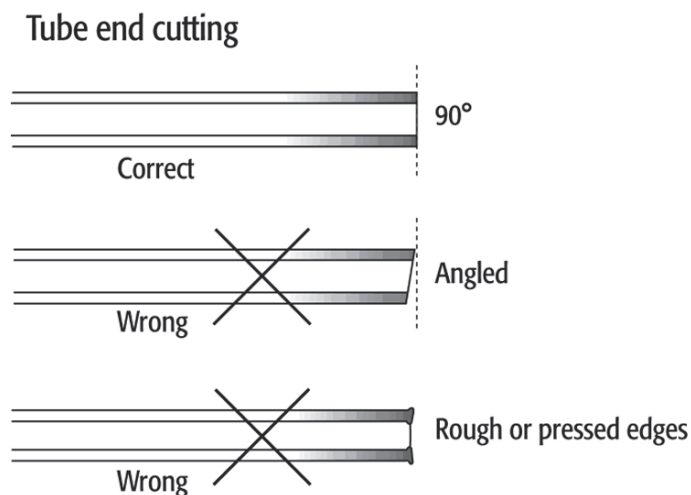
All tubes have to be cut absolutely square (Figure 59). Any deviation from square causes a dead volume, which yields carryover and other chromatographic effects. This rule is valid for stainless steel, PEEK, or Polymer tubing. The material is not decisive, but using the correct tool for each material when cutting a tube is.

Stainless steel tubes are often cut by using pliers. Often, an egg-shaped profile results, which no longer seals and causes dead volume. Dedicated pliers for HPLC tubing are available on the market; nevertheless, for best results, use precut tubing which is cut smooth and clean, and is passivated.

Because polymer tubes, such as PEEK, Teflon, PFA, and so on are soft, you might choose to cut them with any handy tool. Commercially available tubing cutters from many manufacturers are available. If the blade does not provide a clean, right-angle cut, however, use a different technique.

The most reliable and common way to cut tubing of any material is with a cutter that has an adjustable blade. Carry out the initial turns, readjust the blade, turn once more, and adjust again until approximately half of the tubing wall is cut. Hold the tube on the two sides of the cut with flat-nose pliers and twist the tube until it breaks. You can use the same procedure for polymer tubing by replacing flat-nose pliers with tweezers with a flat tip.

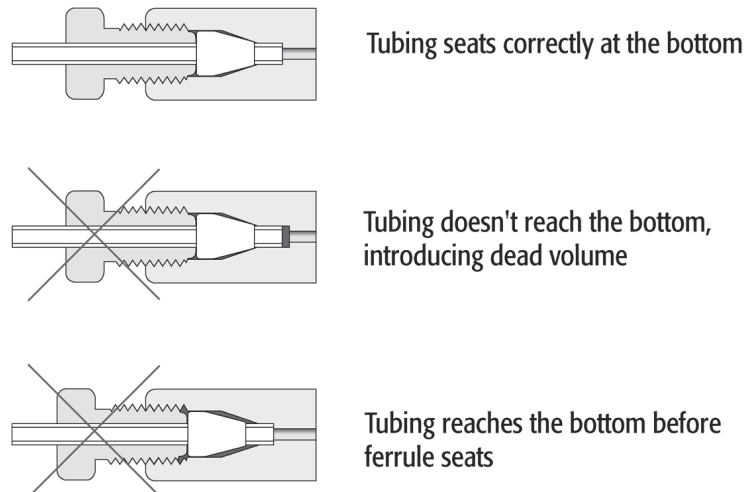
**Figure 59.** Illustration of various tubing cuts



### Wrong Nuts or Ferrules Used

Do not use a nut or ferrule from a vendor other than those specified for the product. [Figure 60](#) illustrates the resulting dead volume when you use an incorrect nut or ferrule type for connecting.

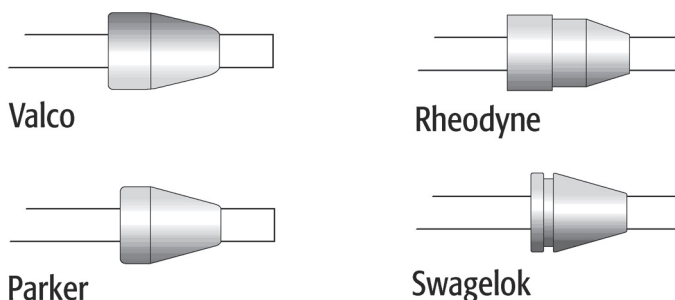
**Figure 60.** Dead volume created due to wrong ferrule type



[Figure 61](#) shows the various forms of the ferrules from different manufacturers. Although some of them are close in shape to what you need, they are not close enough to exchange without risk.



**Figure 61.** Ferrule types from various vendors



### Avoiding Trapped Air Bubbles

It is common practice to make a connection only if the ports are wetted. The liquid helps to prevent trapped air bubbles, which are often tedious to remove or work out of a system.

### Solvent Delivery Pumps Not Priming

The solvent delivery pumps of the Pump Module can fail to prime at initial operation.

You must wet the pump at the solvent inlet port or, preferably, prime the tubing from the Solvent reservoir to the pump and make the connection.

Dry running the pump for a short time will not harm it.

### Solvent Delivery Pump Not Functioning

The pump flow rate depends greatly on the viscosity of the selected wash solvent and the given backpressure of the entire system. Consider the valve bore size (standard Cheminert valve 0.25 mm), as well as the tubing and injection loop internal diameter.

If in doubt about the delivered volume of a pump, run the following test (Test 1). If the result of the test is negative, check the flow from the pump directly (Test 2).

#### ❖ (Test 1) To test the pump flow delivered throughout the entire system

1. Prepare a graduated cylinder with a volume of 10 or 20 mL.
2. Test a solvent, for example, water.
3. Select this path from the local terminal workstation.

**Menu > Utilities > Wash Station > *selected wash station***

The delivery pump and the DLW Actuator are activated, and you can collect the delivered flow at the DLW syringe needle.

Expected result: Collection of approximately 8 mL

(Water, DLW syringe needle, gauge 22)

❖ **(Test 2) To test the pump flow delivered at the pump outlet**

1. Disconnect the outlet tube from the pump module.
2. Connect PEEK tubing with the following dimension to the outlet of the pump module:  
85 mm (3.35 in.) ID, 0.25 mm (0.010 in.) ID.
3. If necessary, replace the solvent in the reservoir bottle with water.
4. Activate the pump by selecting this path:

**Menu > F3-Setup > Objects > Events > Pwr-Out1 or Pwr-Out2**

5. Collect the pumped water in a graduated cylinder.

Expected Result: Within 30 seconds, pump delivery of approximately 10 mL

This short test can indicate whether the system is performing as expected. If you observe no or very little flow, systematically check using a process of elimination. For more details, see [Table 29](#).

## Standard Chromatographic Tests

This section describes standard chromatographic tests for carryover and repeatability. These tests provide basic information for both users and technical support.

The tests described are examples only; vary them accordingly if the configuration of the HPLC system changes.

The first test uses PAL standard valves with a bore size of 0.25 mm to provide a flow rate range of 10–500  $\mu\text{L}/\text{min}$ .

Both tests, carryover and repeatability, test conditions for UV and LC/MS detection.

Details are listed in the next section. Note that the given levels to reach for carryover and repeatability do not reflect the specification value. These described tests should give you confirmation if the system is working to expectation using reasonable effort and time. The tests are more comparable with an OQ Test (Operational Qualification Test).

**Note** The levels provided in these tests do not reflect the specification value for careover or repeatability. The test results will confirm if the system is working properly with reasonable effort and within the expected time frame. The tests are comparable to Operational Qualification (OQ) tests.

## Carryover Test with UV-Detection

Table 31 lists the specifications for carryover tests with UV-detection.

**Table 31.** Specifications for carryover tests with UV-detection

Component	Specification
Test Sample	Chlorhexidine, 0.6 mg/mL dissolved in Water:acetonitrile = 90:10 + 0.1%TFA
Blank Solution	Water:acetonitrile + 0.1%TFA
HPLC System	<ul style="list-style-type: none"> <li>Binary solvent delivery pump</li> <li>UV-Detector, wavelength: 257 nm</li> <li>HPLC column: C18 (Halo), 2.1 x 50 mm spherical 2.7 <math>\mu</math>m</li> <li>Column temperature: Ambient</li> <li>Mobile Phase:               <ul style="list-style-type: none"> <li>A: Water + 0.1%TFA</li> <li>B: Acetonitrile + 0.1% TFA</li> </ul> </li> <li>Gradient:               <ul style="list-style-type: none"> <li>10% A hold for 0.5 min, to 90% B in 7.5 min hold for 1 min</li> </ul> </li> </ul>
Wash Solvents	<ul style="list-style-type: none"> <li>Wash1: Water + 0.1% TFA</li> <li>Wash2: Acetonitrile + 0.1% TFA</li> </ul>
Loop Size	2 $\mu$ L; Loop ID 0.13 mm
Injection Valve	VICI/Valco (CH-106678), bore size 0.25 mm
Cycle Composer Macro or Cycle (ICC-CE)	DLW Standard

### Cycle Composer Method Parameter Settings

Table 32 lists the Cycle Composer Method parameter settings for carryover tests with UV-detection.

**Table 32.** Carryover tests with UV-detection (Sheet 1 of 2)

Parameter	Value
Syringe	100 $\mu$ L DLW
Sample Volume	7 $\mu$ L
Airgap Volume	3 $\mu$ L
Front Volume	5 $\mu$ L
Rear Volume	5 $\mu$ L

**Table 32.** Carryover tests with UV-detection (Sheet 2 of 2)

Parameter	Value
Needle Gap Valve Clean	3 mm
Fill Speed	5 $\mu$ L/s
Fill Strokes	0
Stator Wash	0
Pullup Del	500 ms
Inject to	Specified Inj. Valve
Inject Speed	5 $\mu$ L/s
Pre Inj Del	500 ms
Pst Inj Del	500 ms
Valve Clean Time Solvent 2	2 s
Post Clean Time Solvent 2	2 s
Valve Clean Time Solvent 1	3 s
Post Clean Time solvent 1	2 s

## Repeatability Test with UV-Detection

Table 33 lists the specifications for repeatability tests with UV-detection.

**Table 33.** Repeatability tests with UV-detection (Sheet 1 of 2)

Component	Specification
Test Sample	<ul style="list-style-type: none"> <li>• Loop overflow: Caffeine, 520 mg/L</li> <li>• Partial loop: Caffeine, 52 mg/L dissolved in water:acetonitrile = 90:10 + 0.1%TFA</li> </ul>
HPLC System	<ul style="list-style-type: none"> <li>• Binary solvent delivery pump</li> <li>• UV-Detector, wavelength: 273 nm</li> <li>• HPLC column: Hypersil Gold C18, 2.1 <math>\times</math> 50 mm spherical 1.9 <math>\mu</math>m</li> <li>• Column temperature: Ambient</li> <li>• Mobile phase Water:acetonitrile = 90:10 +0.1% TFA</li> <li>• Gradient: Isocratic for 2 min</li> </ul>
Wash Solvents	<ul style="list-style-type: none"> <li>• Wash1: Water + 0.1% TFA</li> <li>• Wash2: Acetonitrile + 0.1% TFA</li> </ul>

**Table 33.** Repeatability tests with UV-detection (Sheet 2 of 2)

Component	Specification
Loop Size	<ul style="list-style-type: none"> <li>• 2 µL (Loop overfill, 7 µL)</li> <li>• Loop ID: 0.13 mm</li> <li>• 20 µL (partial loop, 10 µL)</li> <li>• Loop ID 0.25 mm</li> </ul>
Injection Valve	VICI/Valco C72VX-1696D-CTC, bore size 0.25 mm
Cycle Composer Macro or Cycle (ICC-CE)	DLW Standard

#### Cycle Composer Method Parameter Settings

Table 34 lists the Cycle Composer Method parameter settings for repeatability tests with UV-detection.

**Table 34.** Repeatability tests with UV-detection

Parameter	Value
Syringe	100 µL DLW
Sample Volume	7/10 µL
Airgap Volume	3 µL
Front Volume	5 µL
Rear Volume	5 µL
Needle Gap Valve Clean	3 mm
Fill Speed	5 µL
Fill Strokes	0
Stator Wash	0
Pullup Del	500 ms
Inject to	Specified Inj. Valve
Inject Speed	5 µL/S
Pre Inj Del	500 ms
Pst Inj Del	500 ms
Valve Clean Time Solvent 2	2 s
Post Clean Time Solvent 2	2 s
Valve Clean Time Solvent 1	3 s
Post Clean Time Solvent 1	2 s



## Replaceable Parts

This chapter describes the replaceable parts for the Accela Open Autosampler and the Dynamic Load and Wash (DLW) option.

### Contents

- [DLW Pump Module](#)
- [DLW Wash Station](#)
- [DLW Syringe](#)
- [DLW Syringe Holder Assembly](#)
- [DLW Option Spare Parts Ordering Information](#)
- [Control-xt Board](#)
- [Injection Valve and Valve Rotor](#)
- [Injection Unit](#)



### CAUTION

- Before attempting any type of maintenance, always disconnect the power cords from the power supply or from the various power supplies if optional devices are installed.
- To prevent possible shock, do not touch the capacitors. Capacitors inside the instrument can remain charged, even if the instrument is turned off.
- To avoid damaging electrical parts, do not disconnect an electrical assembly while power is supplied to the PAL System. Once the power is turned off, wait approximately 30 seconds before you disconnect an assembly.

This section provides short descriptions of parts that can be replaced in the field.

## DLW Pump Module

You can exchange the following parts individually from the DLW pump module:

- Single-pump module (incl. electrical connection)
- Solvent tubing kit
- Electrical cable (from pump module to PAL System)
- Pump control PCB
- DLW tubing kit

## DLW Wash Station

You can exchange the following parts individually from the DLW wash station module.

- Wash station block
- Wash station inserts
- Waste tube
- Waste tube (valve-to-wash station)
- Dummy plugs

**Note** The inserts and individual parts of the wash station block will be available in an update for this system part.

## DLW Syringe

You can exchange the following parts individually from the DLW syringe assembly:

- DLW syringe 100  $\mu$ L
- DLW syringe plunger (for syringe 100  $\mu$ L)
- DLW syringe plunger holder
- DLW syringe retaining nut
- DLW syringe replaceable needles



## DLW Syringe Holder Assembly

You can exchange the following parts individually from the DLW syringe holder assembly:

- DLW syringe holder assembly (complete assembly)
- DLW needle adapter block
- DLW needle holder assembly
- Holding loop
- DLW manifold

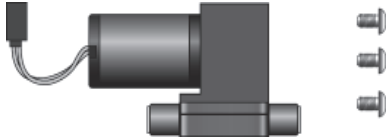
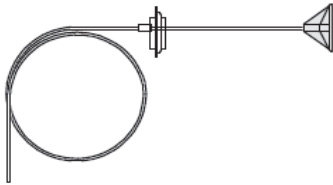
(Entire unit assembled, consisting of DLW solvent selector and DLW syringe adapter with Perfluor O-ring installed)

- DLW flow diverter
- DLW actuator/solenoid
- DLW actuator control PCB (incl. cable)
- DLW syringe holder fixation screw

## DLW Option Spare Parts Ordering Information

Table 35 contains spare parts ordering information for the DLW option.

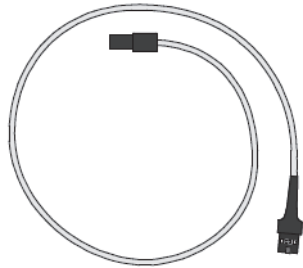

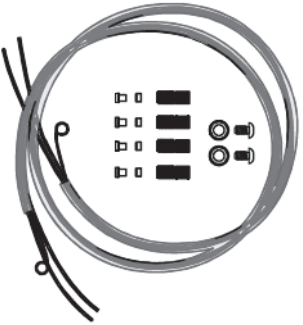
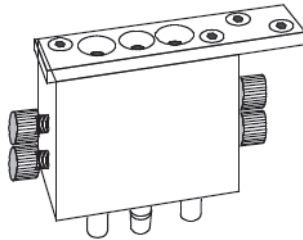
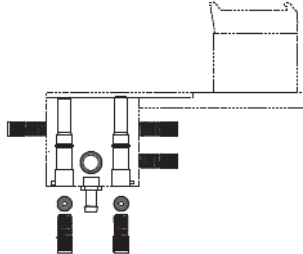
**Table 35.** DLW option spare parts ordering information (Sheet 1 of 5)

Part name/number	Description	Illustration
<b>DLW Pump Module</b>		
00950-01-00310	PAL DLW Pump Kit (single module), incl. electrical connections, 24 V	
00950-01-00311	Aspiration Tube Kit, consisting of: <ul style="list-style-type: none"> <li>• 1 pc. Tube PFA</li> <li>• 1 pc. PEEK solvent filter 10 µm</li> <li>• 1 pc. Plug with feed through hole</li> </ul>	




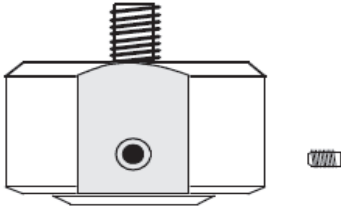
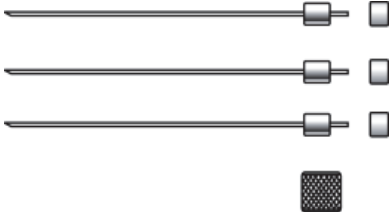
## 7 Replaceable Parts

DLW Option Spare Parts Ordering Information

**Table 35.** DLW option spare parts ordering information (Sheet 2 of 5)

Part name/number	Description	Illustration
00950-01-00312	Cable for PAL-xt wash station round connector, 130 cm long	
00950-01-00313	PAL DLW Pump Control Module PCB Kit	
00950-01-00314	PAL DLW Wash Station Tubing Kit, consisting of: 2 pcs. tubing, wire supported and connections	
<b>DLW Wash Station</b>		
00950-01-00315	PAL DLW Wash Station Block, complete assembly (incl. centering plate)	
00950-01-00334	PAL DLW Wash Station Insert Kit, consisting of: <ul style="list-style-type: none"><li>• 2 pcs. Inserts</li><li>• 1 pc. Waste Connector</li><li>• 5 pcs. Dummy Plugs</li></ul>	

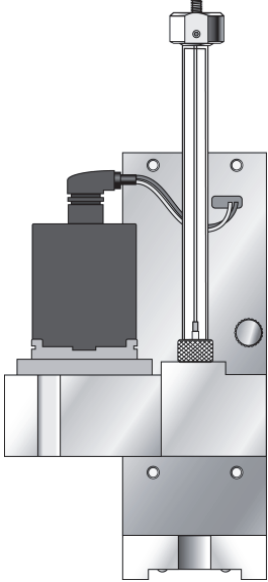
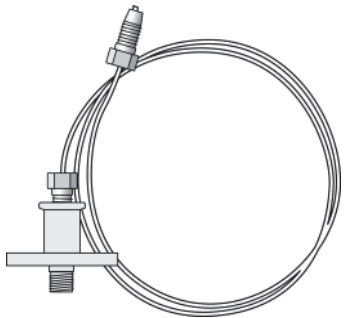
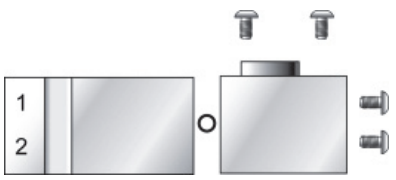

**Table 35.** DLW option spare parts ordering information (Sheet 3 of 5)

Part name/number	Description	Illustration
00950-01-00335	Tube waste for wash station, length 2 m	
<b>DLW Syringe</b>		
00950-01-00325	Syringe for PAL DLW Option, 100 µL gastight; Ø7.7 mm; Scale Length 60 mm; Thread ¼-28 UNF. Removable needle, not included	
00950-01-00326	Replacement plunger for syringe SYRC DLW-R.; pkg of 10	
00950-01-00327	PAL DLW Plunger Holder	
00950-01-00328	Needle Kit for PAL DLW Option, gauge 22 PST 3, / 51 mm (3 pcs. per pack); incl. needle retaining nut	

## 7 Replaceable Parts

DLW Option Spare Parts Ordering Information

**Table 35.** DLW option spare parts ordering information (Sheet 4 of 5)

Part name/number	Description	Illustration
<b>DLW Syringe Holder assembly</b>		
00950-01-00316	PAL DLW Syringe Holder assembly; complete assembly with 100 $\mu$ L syringe, without syringe needle	
00950-01-00317	Kit PAL DLW Holding Loop with needle adapter mounted	
00950-01-00318	PAL DLW Manifold Kit, complete, assembled	
00950-01-00319	PAL DLW Flow Diverter, pkg of 1	

**Table 35.** DLW option spare parts ordering information (Sheet 5 of 5)

Part name/number	Description	Illustration
00950-01-00320	PAL DLW Actuator/Solenoid Kit, incl. screws, without cable	
00950-01-00321	PAL DLW Actuator Control PCB Kit, incl. cable and screw set	
00950-01-00322	Syringe Holder Fixation Screw, pkg of 1	
00950-01-00323	PAL DLW Needle Length Guide tool, pkg of 1	
00950-01-00324	PAL DLW Needle Holder Kit	

**Note** Individual parts or the total composition of spare part kits can change without notice.

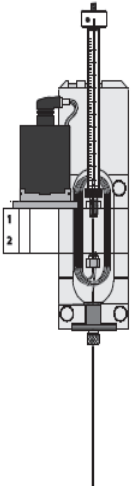
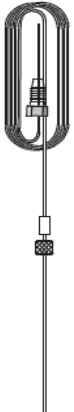
## 7 Replaceable Parts

DLW-2 Option Spare Parts Ordering Information

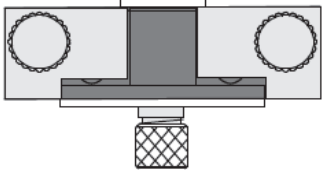
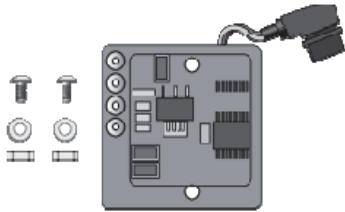
# DLW-2 Option Spare Parts Ordering Information

Table 36 contains spare parts ordering information for the DLW-2 option. The complete list of the DLW-2 spare parts include all DLW spare parts and the parts in Table 36.

**Table 36.** DLW-2 option spare parts ordering information (Sheet 1 of 2)

Part name/number	Description	Illustration
<b>DLW Pump Module</b>		
00950-01-00359	PAL DLW-2 Syringe Holder assembly; complete assembly with holding loop and 100 $\mu$ L syringe installed	
00950-01-00360	Kit PAL DLW-2 holding loop with need and adapter mounted	

**Table 36.** DLW-2 option spare parts ordering information (Sheet 2 of 2)

Part name/number	Description	Illustration
00950-01-00361	PAL DLW-2 Needle Holder Kit	
00950-01-00362	PAL DLW-2 Actuator Control PCB Kit, incl. cable and screw set	

## Control-xt Board

**PN: 00950-01-00329**

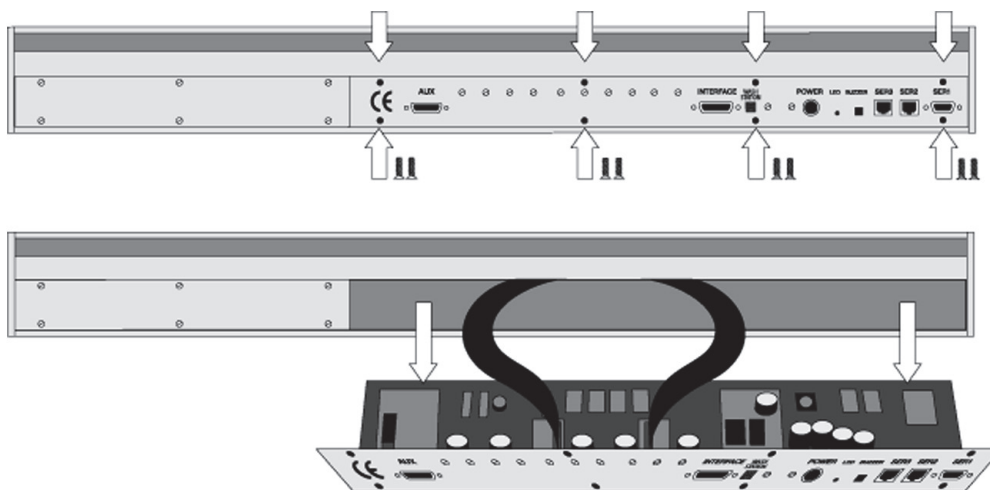
### ❖ To install the control-xt board

**Note** The Control-xt board replaces the control board (PN CH-952589). This board is required for the Accela Open Autosampler system and operates only at PAL Firmware level 4.1.x or later.

1. Follow the sequence shown in [Figure 62](#) to detach and release the Control-xt board from its position inside the *x* axis. As shown, remove the dummy cover first.
2. Carefully pull the two ribbon cable connectors upwards to free them from the board.
3. Install the replacement board in the reverse order.

**Tip** After reconnecting the two ribbon cables to the board, ensure that the cables are folded flat and can slide inside the *x* axis without damage.

**Figure 62.** Replacing Control-xt board



## Injection Valve and Valve Rotor

This section describes the commonly used injection valves with 6-ports.

### Cleaning or Replacing Injection Valves

This section contains information about cleaning or replacing the injection valve for the VICI/Valco Cheminert.

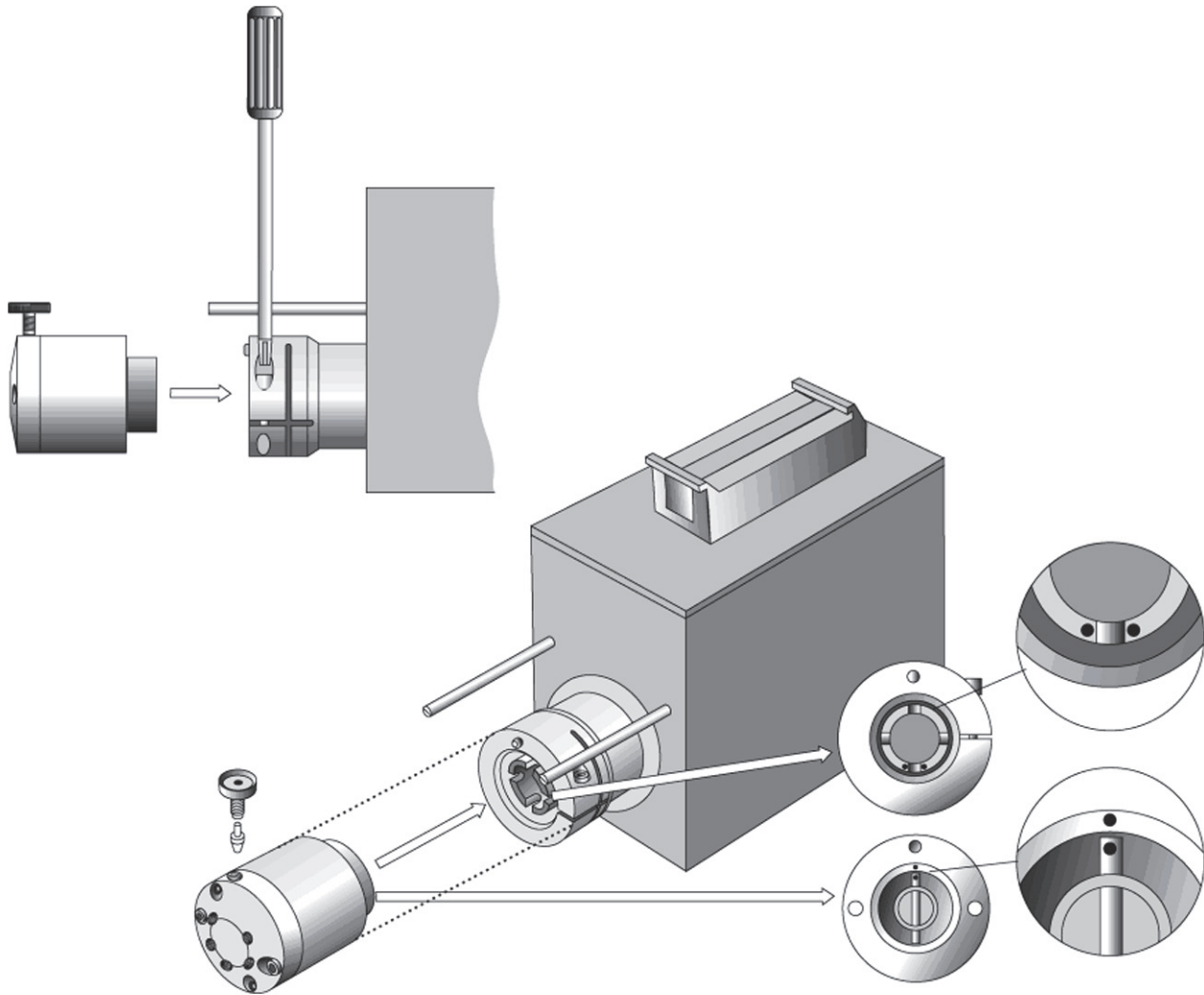
#### Cheminert

❖ **To clean or replace the Cheminert injection valve (Figure 63)**

1. Remove the sample loop, needle guide fitting, and connection tubes from the valve body.
2. Unscrew the Torx screw at the valve adapter, which holds the valve to the valve drive.
3. Check that the valve guide pin remains at its position and observe the guide marks as shown in [Figure 46](#) on [page 50](#).
4. Install the replacement valve in the reverse order.



**Figure 63.** Replacing the VICI/Valco Cheminert injection valve



## Cleaning or Replacing Valve Rotors

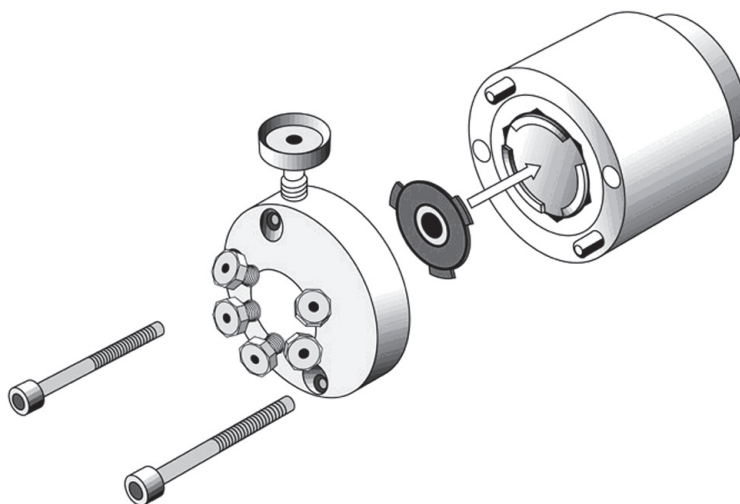
This section contains information about cleaning or replacing valve rotors for the VICI/Valco Cheminert.

### Cheminert Valve Rotor

❖ **To clean or replace the valve rotors for VICI/Valco Cheminert (Figure 64)**

1. Do not remove the valve from the valve drive. For improved access, you may remove all tubing or the loop.
2. Unscrew the two hex screws, which hold the valve stator to the valve body.
3. Carefully lift out the rotor disk with a protected tool (with tips protected to avoid scratching the surface).
4. Insert the new rotor. The size-coded positioning tabs prevent you from inserting the rotor upside down.
5. Tighten the hex screws carefully and evenly to reach a parallel seat of valve stator and body.

**Figure 64.** Rotor replacement of Cheminert valve



## Injection Unit

For details on the parts of the injection unit, see “Installing the Injection Unit” on page 18.

## Reference Information

This appendix contains the following reference information.

### Contents

- [Glossary](#)
- [Naming Convention](#)
- [Special Functions](#)
- [PAL Firmware Overview](#)
- [PAL Firmware Error Codes](#)
- [External Connectors for Accela Open Autosampler](#)
- [System Limitations](#)
- [Interfacing the Open Autosampler to Other Devices](#)

## Glossary

Table 37 provides a list of glossary terms related to the Accela Open Autosampler.

**Table 37.** Accela Open Autosampler glossary (Sheet 1 of 2)

Terms	Description
Cycle	The specific operations necessary to process one sample. The cycle operations are repeated for each sample within a job. Cycles are designed for specific applications.
Job	The information needed by the Accela Open Autosampler to process multiple samples by the same processing steps. The elements of a job are a method and a tray that define the location of the samples to be processed. For identification, jobs are automatically numbered 01–99 and then restart with 01 when they are added to the Job Queue.
Job queue	A list of sample processing jobs. Jobs are executed in the order displayed on the JOB QUEUE menu screens. You can add new jobs to the queue while samples are being processed.
Method	How the samples are processed. The elements of a method are a cycle, a syringe, and a parameter list. Methods have names with up to eight characters and can be edited, copied, and deleted.
Method parameters	Associated with the cycle operations. User-assigned parameter values define how a processing operation is performed.  A zero parameter value will disable a cycle operation. Cycle parameters are application-specific.
Module	PAL hardware module, either part of a standard PAL configuration (for example, Combi PAL, HTS PAL) or an optional addition (for example, cooled stack, MALDI tool, dilutor).  The term <i>Module</i> is intentionally used to differentiate from object, which is reserved for the PAL firmware object.
Objects	Data structures describing the properties of physical modules. Certain modules (for example, a stack) require several objects.
Object class	Each object belongs to an object class (for example, syringes, trays, injectors).  The object class defines the Items of an object.
Object Item	An object contains several Items, which can be numerical values with a physical unit (for example, $x$ -, $y$ -, $z$ -position, penetration, syringe scale length, syringe volume) or references to other objects.  Note that the term, <i>Parameter</i> , is reserved for ATOM Parameter.  (PAL firmware commands are to be used for a PAL cycle or macro).

**Table 37.** Accela Open Autosampler glossary (Sheet 2 of 2)

<b>Terms</b>	<b>Description</b>
PAL Object list	<p>If you add a PAL module (hardware) to an instrument, you must load several objects into the firmware. These objects are collected in an object list and stored in a file with the extension .pol.</p> <p>Object lists are delivered together with object manager software and are grouped into folders for the different kind of modules (for example, syringes, tray holders, valve drives). The name of an object list starts with the module part number with variants added (for example, first or second stack). The name of the root folder includes the revision which depends on the firmware version (for example, object lists Rev. K for firmware 2.x and 3.x).</p>
PAL Object Manager	Software to load a PAL object list to an instrument if a module (hardware module) has been added to the PAL System. In a special mode, you can also use the Object Manager to create and maintain object lists.
Stack	<p>A particular type of tray holder that is designed to hold microplates. A six-drawer stack holds 12 standard microplates, two in each drawer.</p> <p>A three-drawer stack holds six deep-well microplates, two in each drawer.</p>
Tray	Holds multiple samples. Trays are defined by designating the tray type (see below) and the tray holder. Tray names are used to identify the sample source within an autosampler job.
Tray holder	Holds one or more trays. Each tray holder has a reference position ( $x$ , $y$ , $z$ coordinates) that defines its location.
Tray type	The pattern and sampling sequence of sample locations within a tray.

## Naming Convention

This section recommends the standard naming convention for the Accela Open Autosampler. Following this convention will allow the PAL setup to be preconfigured for certain applications. This will simplify software backups and application development, and will improve technical support and training.

**Table 38.** Naming conventions

Tray type	Tray description
VT200	Vial tray, 200 positions (10 × 20)  For 7 mm micro vials 1 mL
VT98	Vial trays, 98 positions (7 × 14)  For 12 mm vials 2 mL
VT78	Vial tray, 98 positions (6 × 13)  For 7 mm micro-vials, 1 mL (opposite side of 98 positions tray)
VT54	Vial tray, 54 positions (6 × 9)  For 12 mm vials, 2 mL
VT21	Vial tray, 21 positions (7 × 4)  For 12 mm vials, 2 mL
VT32-10	Vial tray, 32 positions (4 × 8)  For 23 mm headspace vials, 10 mL
VT32-20	Vial tray, 32 positions (4 × 8)  For 23 mm headspace vials, 20 mL
VT15-10	Vial tray, 15 positions (3 × 5)  For 23 mm vials, 10 mL
MT96	Standard 96-position shallow microplate
DW96	Deep well 96-position microplate
MT384	High density 384-position shallow microplate

## Special Functions

This section describes special functions that you can activate by applying the Extended User mode.

The purpose of having two software access levels is to display discrete items and sections of the firmware at each level. Settings that you set rarely—perhaps at the time of the Accela Open Autosampler system installation—are hidden at level 1 (User Level) and revealed at level 2 (Extended User Level). This also protects the Accela Open Autosampler in group use. Nevertheless, you should become familiar with the important points as described.

### Accessing Extended User Mode

Access the Extended User mode by the following path:



You can navigate to the Extended User mode as follows:

1. Choose **Menu > F3-Setup**.
2. Activate Setup by pressing the center of the selector knob.

For the purpose of the PAL user manual or technical notes, use the following only:

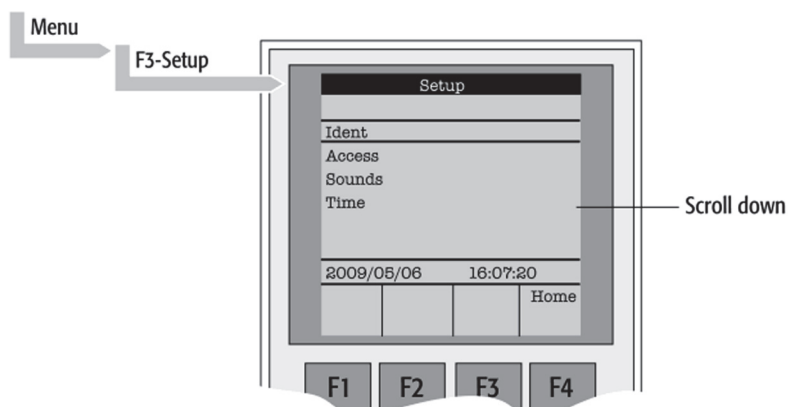
**Menu > F3-Setup.**

### Section F3-Setup

This Setup section at the Extended User Level provides various options to access other classes that are not visible at the User Level. You might already be familiar with some of the classes but others might be new. See the details in [Table 39](#).

To open this level, use the following path ([Figure 65](#)).

**Figure 65.** Selecting Setup in Extended User mode



**Table 39.** Extended User Level options

Object	Description
Ident	The Identification of the system provides the opportunity to assign users, site, and system names. You can enter the serial number of the Accela Open Autosampler at this level. After loading a fresh PAL Firmware backup file, the SNo. is displayed as XXXXXX.
Access	Selectively open or limit access to different users for the Job, Method, or Setup.
Sounds	An adjustable beep sound for specific needs. For safety reasons, keep the default settings, especially the Warn Move, turned on to signal the start of the PAL System to the user.
Time	Sets the PAL internal clock.  If the clock is fast, reset all items, from Year to Seconds and use the mandatory function key F1 (for Set Time).
Communication	Define the communication mode, serial or LAN, and define conditions.  The current Accela Open Autosampler does not support LAN communication.
System	Predefine basic parameters for specific needs.
Service	Activate basic service tests.
Objects	Accessing the PAL Firmware Class Objects in the Extended User mode provides more detailed possibilities for object items. At this level, you can copy or delete an Object by using the specific function keys. This provides quick access to add, for example, a second or third injector.



## System

Select the particular item to change system settings.

PAL Firmware level 4.1.x adds an intermediate step. After selecting System, the new Settings level follows. Activating Settings opens the described items in [Table 40](#).

**Table 40.** System items<sup>a</sup> (Sheet 1 of 3)

Item	Description
Syr Detect	Turns off the syringe detection if the sensor fails or if a special syringe with unknown ID is installed. Turning the syringe identification sensor off displays a dialog box for selecting a syringe identification.
Start Ref	Provides the choice of referencing all axes at the start of the system or suppressing the referencing. Use this function to detect a possible loss of steps.
Stop XY Error	<p>Turns off the automatic recovery of the PAL System after detecting a collision (loss of steps of a stepper motor).</p> <p>Stop XY Error = ON: Automated recovery turned OFF</p> <p>Stop XY Error = OFF: Automated recovery turned ON</p> <p>If recovery is turned off, the PAL System always checks its position before moving to the injection port for injection. This can result in an extra move, crossing over the <i>x-y</i>-axes sensors, if the injection port is at the opposite end of the <i>x</i> axis relative to the sample location, (zero <i>x</i>-axis position in between sample location and injection port).</p> <p>For critical application fields (for example, clinical analyses), turn off the automated recovery; Stop XY Error= ON.</p>
<i>PlgPathCheck</i>	Monitors the plunger travel path to ensure that the plunger tip reaches the syringe zero point in all cases as compared to the original plunger referencing and defining the syringe zero point. If an error is detected, the PAL System stops. This feature is useful if: sample solutions with particles need to be handled, the syringe type is not adequate for the solution composition, or the syringe plunger is blocked or bent.
PlgChnge Pos	Changes a syringe using the Chnge Syr menu function and moves the plunger up to the position specified by this item. For normal use, accept the default values.

**Table 40.** System items<sup>a</sup> (Sheet 2 of 3)

Item	Description
Init Syr at	<p>Initializes the syringe plunger at the Home or Waste position.</p> <p>If samples with a high danger of toxicity are treated, use the Waste position for referencing.</p>
Inj Signal	<p>An autosampler is the master instrument in a chromatographic system. The other components send a Ready signal to enable the autosampler to start. At the moment of injection, the PAL System sends a Start signal to the other components. The Inj Signal item defines the status when the signal is sent. The syringe is filled with the defined sample volume.</p> <p>PlgUp = The Start signal is sent at the moment the plunger starts moving down for injection.</p> <p>PlgDown = The Start signal is sent at the moment the plunger reaches zero point from syringe. (Recommended for large volume injection, but not in combination with Agilent Chemstation.)</p> <p>ValveSw = The Start signal is sent at the moment the injection valve is switched, HPLC technique.</p> <p>PreInj = Start signal sent at a time before injection, as with a sample preparation device started before a GC or LC system. Minus time relative to chromatographic start time.</p>
Vial PrePress	<p>Avoids a vacuum if a relatively large amount of sample is taken out of a small volume vial. If, for example, a volume of 50 µL is specified for injection, the PAL System first takes 50 µL of ambient air and injects this volume into the sealed vial, and causes overpressure in the vial. Then, the sample volume is aspirated without forming a vacuum. This functionality is only valid for liquid sampling, and not valid for the Headspace technique.</p> <p>Starting with PAL Firmware level 4.1.x, this item is also active in combination with Thermo PAL driver.</p>
LC-Inj	<p>The standard cycle for the HPLC technique.</p> <p>For details, refer to Chapter 5 in the <i>Accela Open Autosampler User Guide for LC Devices</i>.</p>
LC-Cut	<p>The cycle for column switching with two valves for the HPLC technique. Control both valve drives through the AUX interfaces.</p> <p>For details, see “<a href="#">PAL Firmware Overview</a>” on <a href="#">page 136</a>.</p>

**Table 40.** System items<sup>a</sup> (Sheet 3 of 3)

<b>Item</b>	<b>Description</b>
GC-Inj	The standard cycle for the GC technique. For details, see “PAL Firmware Overview” on page 136.
GC-InS	The standard cycle for the GC Sandwich technique. For details, see “PAL Firmware Overview” on page 136.
GC-Dual	The cycle for the GC technique that allows the injection of two samples in two different injectors. The start signal is sent after both sample solutions are injected.  For details, see “PAL Firmware Overview” on page 136 and refer to Chapter 5 in the <i>Accela Open Autosampler User Guide for LC Devices</i> .
HS-Inj	The standard cycle for the GC headspace technique.  For details, see “PAL Firmware Overview” on page 136 and refer to Chapter 5 in the <i>Accela Open Autosampler User Guide for LC Devices</i> .
SPME	The standard cycle for the GC SPME technique, solid phase micro extraction, and fiber techniques. For details, see “PAL Firmware Overview” on page 136.
TTS Sotax	These two cycles are dedicated cycles for dissolution applications. The TTS cycle is used for transdermal patches and the Sotax cycle is used for tablet dissolution.

<sup>a</sup> Items displayed in *italics* are available beginning with PAL Firmware level 4.1.x.

**Note** You can select and activate standard injection cycles directly from the Accela Open Autosampler. You must program any deviation from a standard cycle, such as a Cycle Composer Macro or Cycle Editor (ICC-CE) Cycle, for customized requirements.

**Note** In PAL Firmware levels < 4.0 the standard cycles are grouped in the Cycles item.

## Service

Select the particular item to change service settings ([Table 41](#)).

**Table 41.** Service items<sup>a</sup> (Sheet 1 of 2)

Item	Description
Check Motors	<p>Allows every stepper motor to be moved separately. This function is useful to check out a specific motor or to perform an endurance test after a repair.</p> <p>You can select the Start and Stop positions of the test path, as well as the motor Current, Speed, and Acceleration. You can specify a Pause time. A Cycle Limit and Actual Cycle Counter are useful tools for endurance tests.</p> <p>Phase A: Verifies the functionality of the stepper motor. Manually increase the current for tests. Use this item for troubleshooting only.</p> <p>Phase B: Verifies the functionality of the stepper motor. Manually increase the current for tests. Use for troubleshooting only.</p> <p>MicroStep Inx: Moves the motor within a range of up to four full steps by applying the assigned current.</p> <p>Warning: These items are typically used for service. If applied, do not leave the test settings for longer than a few minutes. Leaving the current setting in a critical range can damage the motor.</p>

**Table 41.** Service items<sup>a</sup> (Sheet 2 of 2)

Item	Description
Test Head	<p>Tests the functionality of the PAL injection unit (Head).</p> <p>Two tests are available by using the function keys:</p> <p>F1 Check Plg: Tests the stepper motor of the plunger drive. Performs a sensor check.</p> <p>F2 Check Z: Tests the vial detection sensor and Needle Guide Blocking function.</p> <p>These tests are only required when the Accela Open Autosampler operates erratically. In such cases, contact a Thermo Fisher Scientific representative.</p>
Test Switches	<p>X-Limit = The actual status of the <i>x</i>-axis sensor. Moves the X-Y-Carriage manually across the limit switch to verify sensor activation.</p> <p>Y-Limit = The actual status of the <i>x</i>-axis sensor. Moves the <i>y</i> axis manually to the <i>y</i>-zero point (at <i>x</i> axis), crossing over the limit switch to verify sensor activation.</p> <p>Z-Limit = The actual status of the <i>z</i>-axis sensor. Moves the syringe slider assembly manually up to the zero <i>z</i>-axis point, crossing over the limit switch to verify sensor activation.</p> <p>Aux1 Limit = The actual status of Aux1 limit sensor.</p> <p>Aux2 Limit = The actual status of Aux2 limit sensor.</p> <p>Plg –Pos = The actual status of Plg–Pos.</p> <p>Needle Guide = The actual status of needle guide sensor.</p>

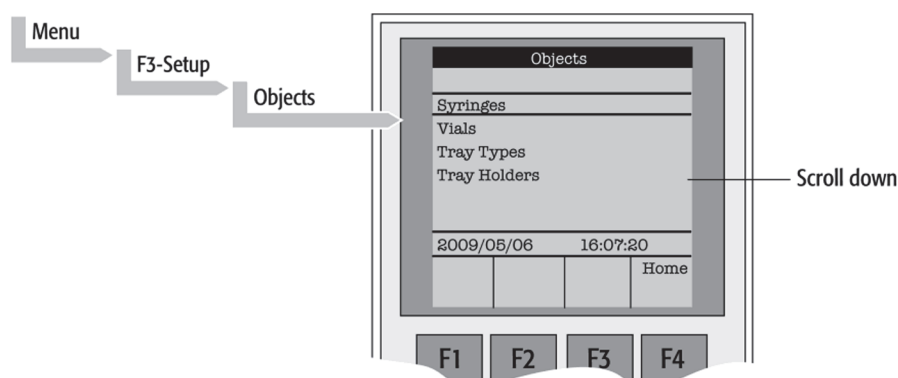
<sup>a</sup> Items displayed in *italics* are available beginning with PAL Firmware level 4.1.x.

## Section F3-Setup/Objects

Use the Objects section opened in the Extended User mode to visualize new classes and known classes that might contain new items accessible only at the Extended User level.

To open this level, use the following path (Figure 66).

**Figure 66.** Selecting Objects in Extended User mode



**Table 42.** Objects in the Extended User mode<sup>a</sup> (Sheet 1 of 4)

Object	Description
Syringes	All loaded syringes are selectable. At the Utilities User Level only the active syringe is visible.
Vials	The Firmware Objects accessed on the PAL Extended User mode have more specific items selectable than at the User Level. Tolerances or ranges such as Maximum Penetration can be defined and limited for the user so that this value cannot be exceeded.
Tray Types	The object class Tray Type is only accessible via the PAL Extended User Level. All loaded Tray Types are visible and selectable.
Tray Holders	The Firmware Objects accessed on the PAL Extended User Level have more specific items selectable than at the User Level. Depending on the tray holder type, an Access and Restore Path or a Heater and Agitator (motor) is assigned.
Trays	All loaded trays are visible and accessible. A tray must be assigned to a tray holder. A TrayTypeGroup identification checks if the specified tray type is allowed in this combination.  The path offset $x$ , $y$ , $z$ is used only if a path is assigned to the specified tray holder. Typically a stack or agitator uses a path.
Wash Stations	The wash station and its positions, Wash1 and Wash2, are visible and selectable.

**Table 42.** Objects in the Extended User mode<sup>a</sup> (Sheet 2 of 4)

<b>Object</b>	<b>Description</b>
Injectors	The Firmware Objects accessed on the PAL Extended User mode have more specific items selectable than at the User Level. Tolerances or ranges such as Maximum Penetration can be defined and limited for the user so that this value cannot be exceeded.
Valves	The valves and valve drives connected and controlled by an AUX interface are listed in this Object class.  The valves controlled by serial communication (serial valves or multiposition valves) are in the PAL Firmware Object class Serial Valves.
Agitators	Items such as Motor Drive, Minimum Speed, and Maximum Speed are selectable for the Agitator Object. You can set the range for the speed. The other items are available directly as Method parameters.
Dilutors	The Firmware Objects accessed in the PAL Extended User mode have more specific items selectable than at the User Level. Tolerances or ranges such as Maximum Fill Speed can be defined and limited for the user so that this value cannot be exceeded.  Items such as Prime Volume, Waste to, Motor Drives, or Scale Length are specific items that remain hidden at the User Level.
Tools	The object class Tools is used to handle another device, such as the MHE-Tool (Multiple Headspace Extraction Tool) or MALDI Tool.
Positions	The various positions used with the PAL System are selectable at this level. Positions such as Home, Change Syringe, or a Path Point, reference the Paths for a specified tray holder directly, such as Agitator or Stack.  Selecting an item (for example, Home) enables the F1 function key and displays Check Pos on the screen. Activating this function provides the possibility of verifying <i>x</i> , <i>y</i> -, <i>z</i> -axes values for this particular position.

**Table 42.** Objects in the Extended User mode<sup>a</sup> (Sheet 3 of 4)

<b>Object</b>	<b>Description</b>
Paths	<p>A Path defines specific movements for the injection unit, which represents a repetitive task assigned to a tray holder. Typical examples are opening and closing a drawer from a Stack or opening the lid of an Agitator.</p> <p>The Path is composed of several steps (points) that are added up in a sequence. Each individual step (point) can be positioned by teaching <i>x</i>, <i>y</i>, <i>z</i> axes. In most cases the return path is in the reverse order of the starting path, for example, AgiOpnL and AgiClsL. If one setting of a point is changed, this change is valid for the return path as well.</p>
Sync Signals	<p>At this level no other functionality, as shown at the User Level, is provided.</p>
Out Signals	<p>At this level no other functionality, as shown at the User Level, is provided.</p>
Events	<p>At this level no other functionality, as shown at the User Level, is provided. Three events are available at this Extended User Level:</p> <p>Pwr-Out1, Pwr-Out2, and FlushVlv. The main purpose of these power-out signals is to activate a solenoid, such as for the wash station or the gas flush valve. You can test the functionality of the signal at this level.</p>
Serial Valves	<p>Serial valves, or multiposition valve drives are powered and controlled by the serial (RS232) control and not through the AUX interface. Starting with firmware level 4.1.x, you can select a mode. The mode refers to the setting of the valve drive, whether the current for the motor is set to High Speed or High Torque.</p>
Out Exp Box	<p>The Out Expansion box provides 8 TTL contacts and 8 relay contacts (24 V contact closure). The optional module (box) is connected and controlled by Interface 2.</p>



**Table 42.** Objects in the Extended User mode<sup>a</sup> (Sheet 4 of 4)

Object	Description
RS232	<p>Use this object to select the serial ports for specific needs. The choices are:</p> <p>Remote, Terminal, Printer, Barcode, ExtDev, VICI-Vlv (for serial valve drives), or None.</p>
<i>Motors</i>	<p>In the class Motors the object Motor-Y is accessible in order to select the items Strategy and Path Speed.</p> <p>Three different settings are selectable for the Y-Motor movement strategy:</p> <ul style="list-style-type: none"> <li>• <b>Rectang</b> = normal movement as used for PAL System. (On firmware level &lt; 4.1.x, this mode was called Auto.)</li> </ul> <p><i>XY-Simul</i> = <i>x</i>, <i>y</i> axes move simultaneously, which allow movement in a diagonal direction. (At firmware level &lt; 4.1.x this mode was called Rectang, but it was not active at lower levels.)</p> <ul style="list-style-type: none"> <li>• <b>Retract</b> = <i>y</i> axis moves from any point first to zero Y (back to the <i>x</i> axis) before the X-movement is started. This mode allows bypassing an object (such as a large GC detector), which might be in the travel path if using the default mode Rectang.</li> <li>• <b>Path Speed</b> = <i>y</i>-axis motor speed in conjunction with a Path, example: Opening and closing a drawer of a Stack. In certain situations you will need to slow down the speed to open and close a drawer. A typical application example is fraction collection, collecting directly into a well plate. If the liquid level is high, the liquid can wash over into the next row if drawer movement is too fast.</li> </ul> <p>Remark: At PAL Firmware level 4.1.x, these items of the Y motor have been made available at the Extended User Level.</p>

<sup>a</sup> Items displayed in *italics* are available beginning with PAL Firmware level 4.1.x.

## Tray Type

You can change Tray Type items by selecting the particular item (Table 43). First select the desired Tray Type, such as **VT32-20**.

**Table 43.** Tray type parameters<sup>a</sup> (Sheet 1 of 3)

Item	Description
Row Length X	Defines the Row length in the X-direction from the center of the first to the center of the last vial position, measured in mm.
Row Length Y	Defines the Row length in the Y-direction from the center of the first to the center of the last vial position, measured in mm.
Col Length X	Defines the Column length in the X-direction from the center of the first to the center of the last vial position, measured in mm.
Col Length Y	Defines the Column length in the Y-direction from the center of the first to the center of the last vial position, measured in mm.
Spl Per Row	Number of samples per Row.
Spl Per Col	Number of samples per Column.
<i>Pattern</i>	<p>The pattern of the Row or Column arrangement has to be defined.</p> <ul style="list-style-type: none"> <li>• Regular = square arrangement (standard)</li> <li>• Staggrd+ = Rows are arranged in a staggered pattern. Offset of second Row is shifted by +50% of hole pattern.</li> <li>• Staggrd- = Rows are arranged in staggered pattern. Offset of second Row is shifted by -50% of hole pattern.</li> </ul>
TrayTypeGroup	Software protection so that not every tray type can be placed on any tray holder. Example: You cannot fit VT32-20 into a Stack.
Plate Thickn	Thickness of tray plate bottom (not the total height of the tray). Measured in mm.
Vial Height	Total height of the vial including cap to seal if applicable. Measured in mm.
	<p>Remark: The sum of the plate thickness and vial height adds up to the total height This is the point where the injection unit expects an object.</p>

**Table 43.** Tray type parameters<sup>a</sup> (Sheet 2 of 3)

Item	Description
Vial Trnsprt	<p>The mode to transport a vial is defined by:</p> <ul style="list-style-type: none"> <li>• <b>None</b> = No transport required. Example liquid sampling.</li> <li>• <b>Magnet</b> = Magnetic Transport. Requires magnetic vial caps. The injection unit moves to the side to slide the vial off when moved away.</li> <li>• <b>Needle</b> = Device is transported with the syringe needle inserted. Example: special vials where a magnetic cap cannot be crimped, for example, Vacutainer.</li> </ul> <p>Remark: This item became active beginning with PAL Firmware level 2.5.x.</p>
<i>ZSlideOffRetr</i>	<p>If magnetic transport is selected, the distance can be defined to move the syringe slider up (Z-direction), lifting up the vial by magnetic force. Measured in mm.</p>
<i>YSlideOffDist</i>	<p>If magnetic transport is selected, the distance can be defined to move the injection unit to the side, Y-direction, to slide the vial off. Measured in mm.</p>
Barcode	<p>If a Barcode Reader is installed, the mode has to be set for <b>AutoFix</b>.</p> <p>None = No Barcode Reader option requested for this tray type.</p>
Z Tolerance	<p>A tolerance window to give a plus/minus range (expressed in mm) where the injection unit must expect an object.</p> <p>If the value of this item is &gt; 0, this is a relative detection mode and the sensor from the injection unit needle guide is active.</p> <p>If the value of this item is set to 0 the sensor of the needle guide is turned off and the syringe slider moves to an absolute value, z-axis position as specified in the object.</p>
Z Retract	<p>The distance to lift up the syringe slider before the y-, z-axes assembly moves across the unit can be specified, expressed in mm.</p>

**Table 43.** Tray type parameters<sup>a</sup> (Sheet 3 of 3)

<b>Item</b>	<b>Description</b>
Max Penetr	<p>The maximum allowed Needle Penetration can be defined as a safety so that the user cannot destroy the needle tip by going too far down.</p> <p>The tray type is prepared for a specific vial size; this value is therefore unique to this combination. The maximum needle penetration must not exceed the length of the syringe needle, considering the loss due to the needle guide, and so on. Typical needle length is 51 mm, and maximum penetration must not exceed 48 mm.</p>
Needle Penetr	This item is identical to that used on the User Level or, if available, as a method parameter of a local cycle. The value defined on these levels is mirrored to the Extended User Level.
Spl1 Offset X	The X offset is defined as the distance from the outer edge of the tray in the X-direction to the center of the vial position 1. Measured in mm.
Spl1 Offset Y	The Y offset is defined as the distance from the outer edge of the tray in Y-direction to the center of the vial position 1. Measured in mm.
Spl1 Offset Z	The Z offset can be selected for special cases where, for example, the syringe needle has to reach a deeper point before penetration into a vial. A Z offset value = 0 considers the tray surface or vial top as the reference point. Measured in mm.

<sup>a</sup> Items displayed in *italics* are available beginning with PAL Firmware level 4.1.x.

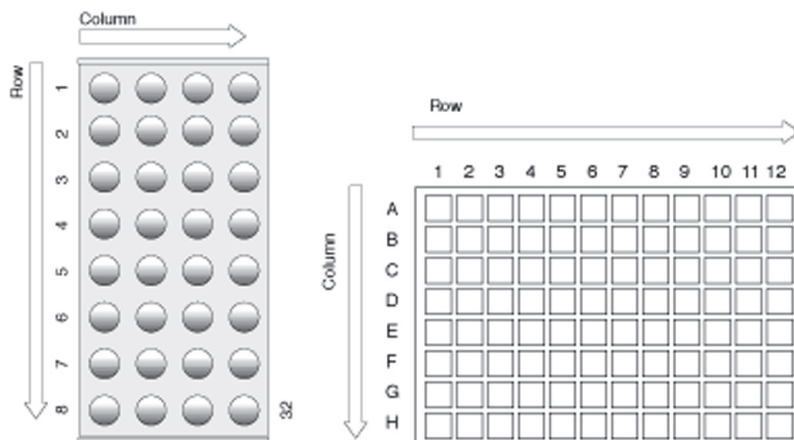
**Note** The dimensions used for Row/Column Length or X-, Y-, Z-Offsets for Sample Number 1, are specific for a particular tray type. If you must make a general adjustment for a tray from a specific vendor—for example, for a Deepwell Plate, make the change in the Tray Type class of Object. For fine tuning, you can use the Offset items from the class Trays.

## Explanation for Various Patterns in Tray Type

- Pattern Type, Regular (Figure 67)

The sample or well positions are equally distributed.

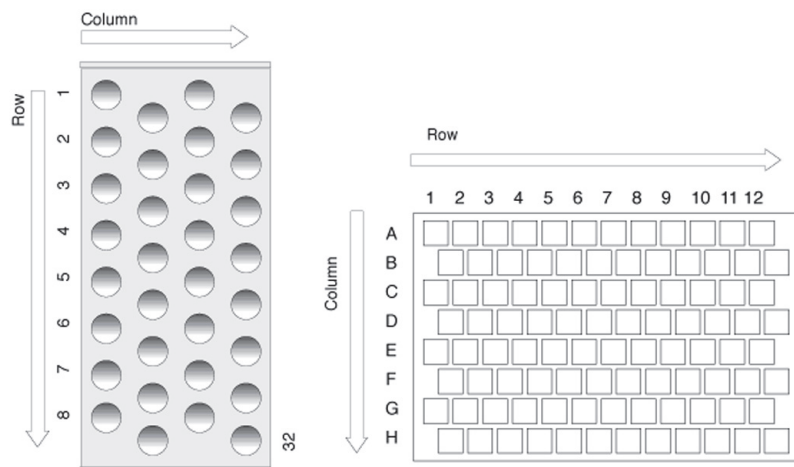
**Figure 67.** Pattern tray type of Regular



- Pattern Type, Staggrd+ (Figure 68)

The sample or well positions are arranged in a staggered pattern. Offset of second Row is shifted by + 50% of hole pattern.

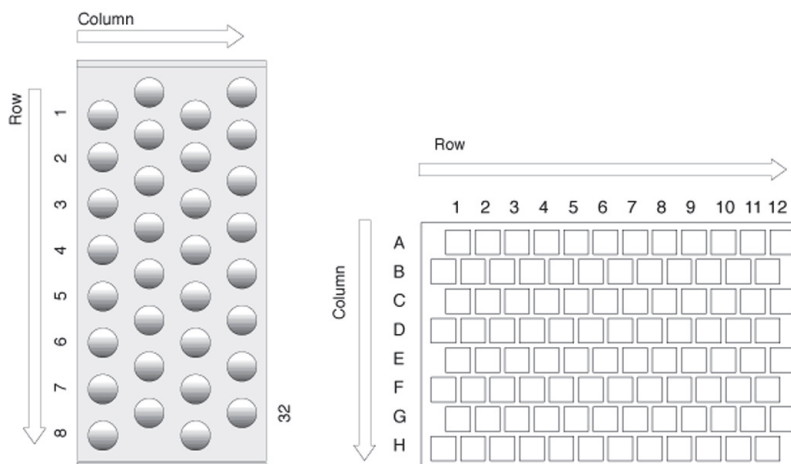
**Figure 68.** Pattern tray type -of Staggered+



- Pattern Type, Staggrd– (Figure 69)

The sample or well positions are arranged in a staggered pattern. Offset of second Row is shifted by –50% of the hole pattern.

**Figure 69.** Pattern tray type of Staggered–



## Trays

You can change the following object Tray Items by selecting the particular item (Table 44). First select the desired tray, such as **Tray1**.

**Table 44.** Tray items<sup>a</sup> (Sheet 1 of 3)

Item	Description
Tray Holder	Specifies the tray holder that the tray is assigned to.
TrayTypeGroup	Displays all tray type groups that are allowed for this particular combination of tray and tray holder. The TrayTypeGroup for the tray type itself is specified for each tray type.
Tray Type	Links the tray type to be used to the tray. This item is usually set at the User Level in Utilities/Trays.
Offset X	<p>The X offset is defined as the distance from the center of the teaching position from the tray holder to the edge of the tray in X-direction. Measured in mm.</p> <p>This item is available at the User Level in the section, Utilities/Tray, and can be used for fine tuning a tray installed in a tray holder. Important for well-plates 384 installed in a Stack.</p> <p>Remark: Position#1 is the only correction point.</p>

**Table 44.** Tray items<sup>a</sup> (Sheet 2 of 3)

<b>Item</b>	<b>Description</b>
Offset Y	<p>The Y offset is defined as the distance from the center of the teaching position from the tray holder to the edge of the tray in Y-direction. Measured in mm.</p> <p>This item is available at the User Level in the section, Utilities/Tray, and can be used for fine tuning a tray installed in a tray holder. Important for well-plates 384 installed in a Stack.</p> <p>Remark: Position#1 is the only correction point.</p>
Offset Z	<p>The Z offset is defined as the distance from the plate surface of the tray holder to the teaching point, the black needle guide that is flush underneath the tray holder plate (which corresponds to the tray holder plate thickness). Measured in mm.</p> <p>This item is available at the User Level in the section, Utilities/Tray, and can be used for fine tuning a tray installed in a tray holder. Important for well-plates 384 installed in a Stack.</p> <p>Remark: Position#1 is the only correction point.</p>
Path Offset X	<p>The Path Offset X can be specific for a tray, adjustment in X-direction. Expressed in mm.</p> <p>Remark: This item is only valid if a path is assigned to the tray holder as specified in this combination.</p>
Path Offset Y	<p>The Path Offset Y can be specific for a tray, adjustment in Y-direction. Expressed in mm.</p> <p>Remark: This item is only valid if a path is assigned to the tray holder as specified in this combination.</p>
Path Offset Z	<p>The Path Offset Z can be specific for a tray, adjustment in Z-direction. Expressed in mm.</p> <p>Remark: This item is only valid if a path is assigned to the tray holder as specified in this combination.</p>
<i>dxRow</i>	Correction for inclination of a tray in X-row direction. For details, see Section 4.2.3.2 above.
<i>dYRow</i>	Correction for inclination of a tray in Y-row direction. For details, see Section 4.2.3.2 above.
<i>dzRow</i>	Correction for inclination of a tray in Z-row direction. For details, see Section F, “Description and Installation,” point 4.2.3.2 above.

**Table 44.** Tray items<sup>a</sup> (Sheet 3 of 3)

Item	Description
<i>dxCol</i>	Correction for inclination of a tray in X-column direction. For details, see Section F, “Description and Installation,” point 4.2.3.2 above.
<i>dyCol</i>	Correction for inclination of a tray in Y-column direction. For details, see Section F, “Description and Installation,” point 4.2.3.2 above.
<i>dzCol</i>	Correction for inclination of a tray in Z-column direction.

<sup>a</sup> Items displayed in *italics* are available beginning with PAL Firmware level 4.1.x.

## Wash Stations

You can change the following object Wash Station items by selecting the particular item (Table 45). First select the desired wash station, such as **Wash1**.

**Table 45.** Wash station items<sup>a</sup> (Sheet 1 of 2)

Item	Description
Position X	<i>x</i> -axis position for the wash station
Position Y	<i>y</i> -axis position for the wash station
Position Z	<i>z</i> -axis position for the wash station
Z Tolerance	<p>A tolerance window for a plus/minus range (expressed in mm) where the injection unit has to expect an object.</p> <p>If the value of this item is &gt; 0, a relative detection mode occurs and the sensor from the injection unit needle guide is active.</p> <p>If the value of this item is set to 0, the sensor of the needle guide is turned off and the syringe slider moves to an absolute value, <i>z</i>-axis position as specified in the object.</p>
Z Retract	Specifies the distance to lift up the syringe slider before the <i>y</i> -, <i>z</i> -axes assembly moves across the unit, expressed in mm.
Max Penetrat	<p>Defines the maximum allowed Needle Penetration as a safety so that the user cannot destroy the needle tip by going too far down.</p> <p>The maximum needle penetration must not exceed the length of the syringe needle, considering the loss due to the needle guide and so on. Typical needle length is 51 mm; maximum penetration must not exceed 48 mm.</p>
Needle Penetr	This item is identical to that at the User Level. The value defined at this level is mirrored in the Extended User Level.



**Table 45.** Wash station items<sup>a</sup> (Sheet 2 of 2)

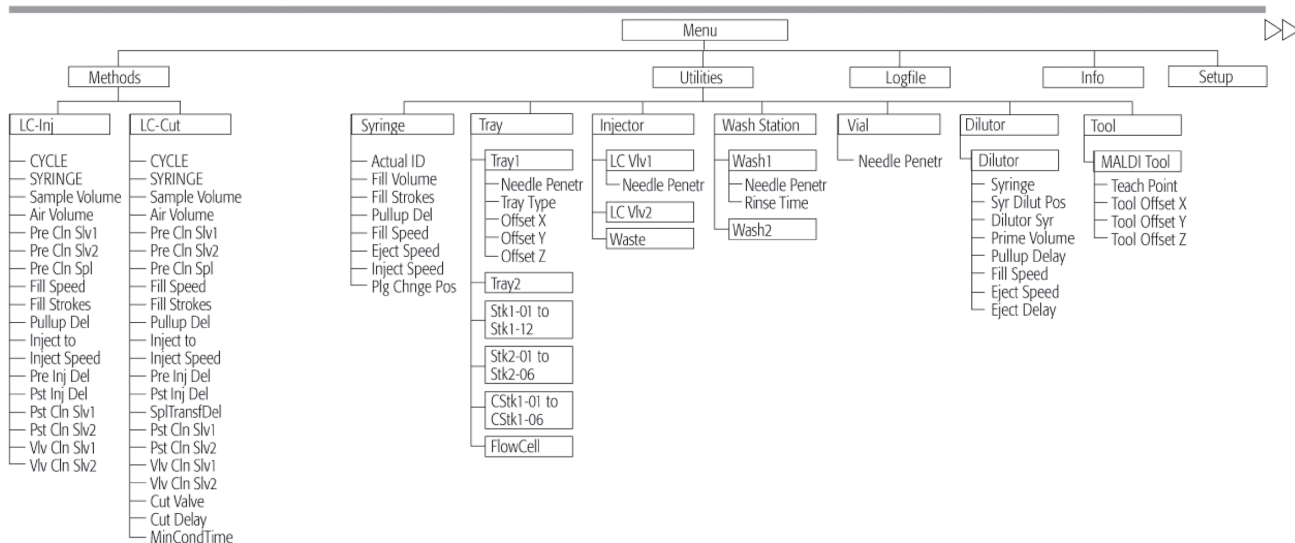
<b>Item</b>	<b>Description</b>
Type	<p>The specific mode for the wash station in use has to be defined:</p> <ul style="list-style-type: none"> <li>• <b>Standard</b> = Standard wash station. Syringe is filled with wash solvent and expelled into Waste.</li> <li>• <b>Pulse</b> = A pulsed signal is sent to activate the device. Used for the fast wash station and active wash station.</li> <li>• <b>Flow</b> = Continuous flow of liquid by gravity. No electrical contact required for a solenoid.</li> <li>• <b>DLW</b> = Dynamic Load and Wash. Item used to activate specific solenoid for DLW Wash Station.</li> </ul>
Flow Control	Specifies the power source to activate a wash station. Uses the Event Power-Out1 and Power-Out2 for the Fast, Active wash station and DLW.
Ndl Cln Path	To assign a path to the wash station.
Waste to	To assign the Waste position to the wash station.
Clean Volume	Defines a percentage of the syringe maximum volume for the cleaning step.
Clean Count	The counter monitors the number of wash cycles. The setting in the method has the same functionality and the value is mirrored in the Extended User Level.
Max Fill Spd	Defines the fill speed for the syringe. If the maximum value of the Fill Speed (10.0 mL/s) is selected, the system takes the injection syringe Fill Speed as defined in the method.
Max Eject Spd	<p>The Eject Speed is the speed used to eject the wash solvent during the wash cycle. This speed is usually higher than the Fill Speed.</p> <p>If the maximum value of the Eject Speed (10.0 mL/s) is selected, the system takes the injection syringe Eject Speed as defined in the object class Syringe.</p>
Rinse Time	<p>The time to rinse the wash port with wash solvent after the wash cycle of the syringe is finished.</p> <p>In the case of a fast/active wash station, the syringe needle is pulled out of the wash port. The wash solvent flows without restriction. As a result, you must consider higher solvent consumption.</p>

<sup>a</sup> Items displayed in *italics* are available beginning with PAL Firmware level 4.1.x.

# PAL Firmware Overview

**Figure 70.** PAL Firmware Overview, Page 1

## Firmware Overview



**Notes:**

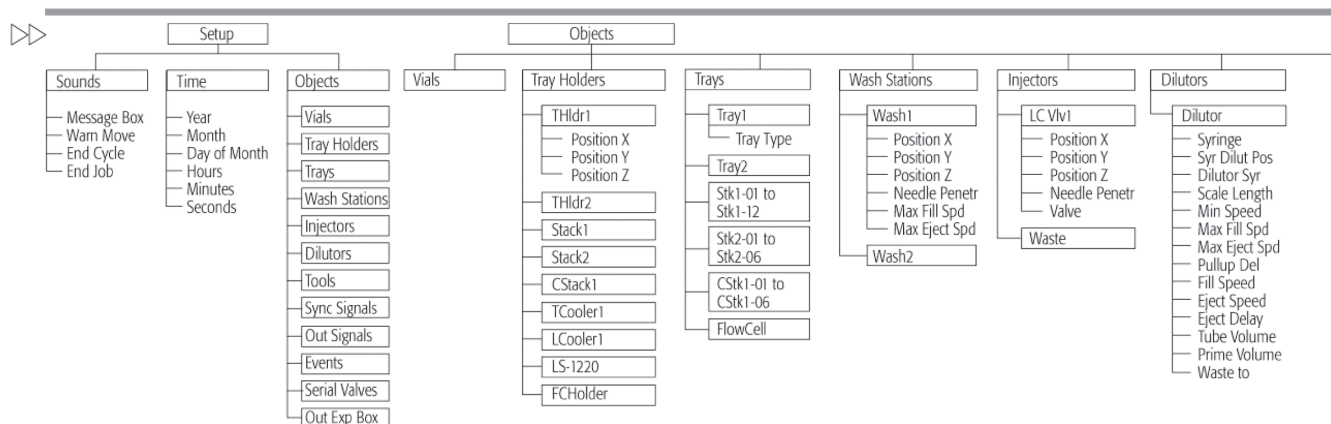
Plunger Stroke and Valve Switch Counter  
Menu/Info/Maintenance

The standard software does not include every Object as shown in the overview.  
The layout depends on the hardware configuration for each individual PAL-System

Revision H / Firmware 4.0.X / June 2009

**Figure 71.** PAL Firmware Overview, Page 2

## Objects



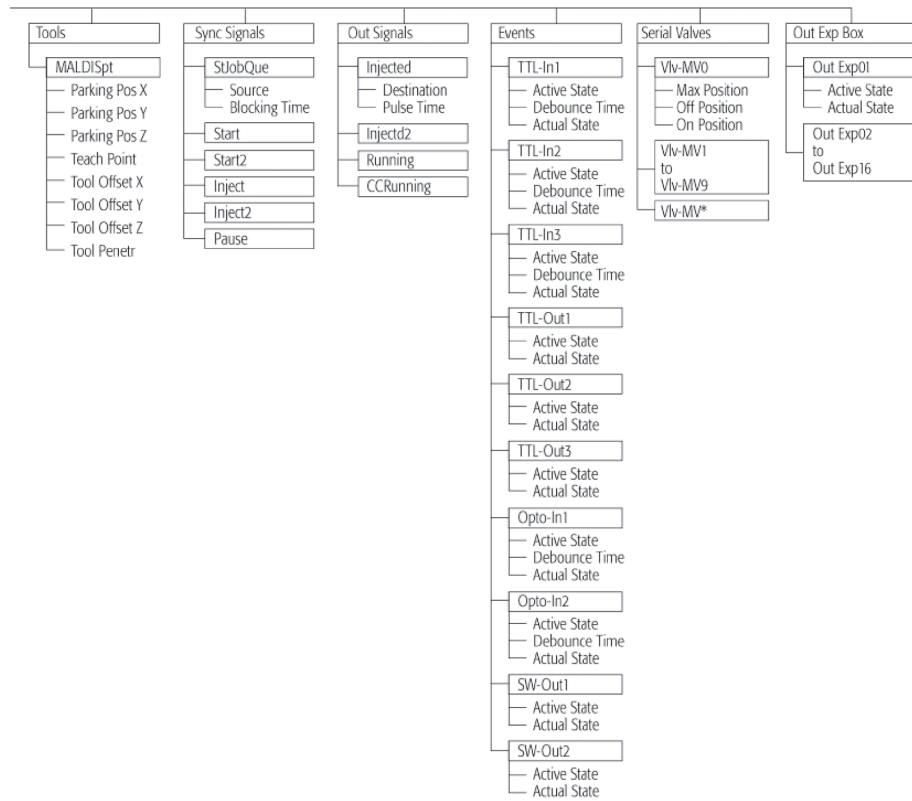
**Note:**  
The standard software does not include every Object as shown in the overview.  
The layout depends on the hardware configuration for each individual PAL-System

Revision H / Firmware 4.0X / June 2009

## A Reference Information

### PAL Firmware Overview

**Figure 72.** PAL Firmware Overview, page 3



PAL is a registered trademark of CTC Analytics AG | Switzerland

**Figure 73.** PAL Firmware Overview, Page 4

Method Cycle	Recommended PAL Method Parameters
<b>LC-Inj</b>	
CYCLE	LC - Inj
SYRINGE	100 µl
Sample Volume	80 µl
Air Volume	0 nl
Pre Cln Slv1	1
Pre Cln Slv2	0
Pre Cln Spl	2
Fill Speed	10 µl/s
Fill Strokes	3
Pullup Del	3-10 s
Inject to	LC Vlv 1
Inject Speed	10 µl/s
Pre Inj Del	500 ms
Pst Inj Del	500 ms
Pst Cln Slv1	2
Pst Cln Slv2	0
Vlv Cln Slv1	2
Vlv Cln Slv2	0

Example for the following conditions:

- Loop Size 20 µL, overfill 3 - 5 times  
Partial Loop filling: Allowed sample volume range 20 - 60% of loop content for loops ≤ 100 µL  
Larger Loops: 20 - 80% of loop content
- Syringe:  
Syr X G100 - 225 - 3: Gauge 225 limits fill speed to max 20 µL/s (5-20 µL/s depends on viscosity of solvent)  
Syr X G100 - 22 - 3: Gauge 22 allows fill speed up to 200 µL/s (depends on viscosity of solvent)
- Eject speed for 100 µL Syringe: 50 to 150 µL/s (Utilities/Syringe)
- Pre - and post Washing:  
Use Solvent 1 and 2 for samples with components of extremely different polarities.  
Samples containing proteins should not contact organic solvents.
- Wash Steps for biological samples:  
1st Wash Cycle: Aqueous solvent  
2nd Wash Cycle: Organic solvent  
  
1st Wash Cycle before next sample:  
Pre-wash with aqueous solvent (eliminate organic solvents in Syringe and Valve)

Method Cycle	Recommended PAL Method Parameters	Remarks
<b>LC-Cut</b>		
CYCLE	LC - Cut	
SYRINGE	100 µl	
Sample Volume	80 µl	
Air Volume	0 nl	
Pre Cln Slv1	1	
Pre Cln Slv2	0	
Pre Cln Spl	2	
Fill Speed	5 µl/s	
Fill Strokes	3	
Pullup Del	3-10 s	
Inject to	LC Vlv 1	
Inject Speed	5 µl/s	
Pre Inj Del	500 ms	
Pst Inj Del	500 ms	
SplTransfDel	Sample Transfer Delay	Time needed to transfer sample from Loop onto 2nd valve (pre - or analytical column)
Pst Cln Slv1	2	
Pst Cln Slv2	0	
Vlv Cln Slv1	2	
Vlv Cln Slv2	0	
Cut Valve	LC Vlv 2	Specify name of 2nd valve, switching valve
Cut Delay		Time needed for clean-up. Switch to backflush sample, to analytical column or detector
MinCondTime		Minimum Conditioning Time: Time needed to condition pre - column on 2nd valve

LC-Cut Cycle controls 2 valves.

- LC Valve 1: Injection valve with loop.
- LC Valve 2: Switching valve.
- Example 1: Pre - column for sample clean-up, backflushing to analytical column. Cut Delay > 0.
- Example 2: 10 - port valve with 2 analytical columns. Loading column 1, condition column 2.  
Valve toggle after Cycle time. Loading column 2, condition column 1. Cut Delay = 0

Part No.: PAL\_FirmOverHrTCr-H

## PAL Firmware Error Codes

Table 46 lists PAL firmware error codes for firmware version 4.1.x.

**Table 46.** PAL firmware error codes (Sheet 1 of 6)

Error code number	Error message
0	Okay = No Error
1	FPGA not present error
2	FPGA configuration error
3	External Event is INPUT only
4	Items list is full
5	Duplicate Object Name found
6	Table String is empty
7	Table String Index is invalid
8	Table String is too long
9	Object list is full
10	Invalid Terminal Firmware
11	Object Name not found
12	Real Time Clock bad Data read
13	Number to String Conversion Error
14	String to Number Conversion Error
15	Object Level too deep
16	Too many Items per Object
17	Item Name not found
18	Axis Position Error
19	Module Trigger not valid
20	Module CRC not valid
21	MenuProc Index out of Range
22	MenuProc Address contains NIL Pointer
23	FuncProc Address contains NIL Pointer
24	Invalid Function Key
25	Invalid MenuTable Index
26	MenuTable Pointer contains NIL
27	Invalid Object Focus in GetItemPtr

**Table 46.** PAL firmware error codes (Sheet 2 of 6)

<b>Error code number</b>	<b>Error message</b>
28	Invalid Item Focus in GetItemPtr
29	Invalid Select String Array Index
30	Unknown Select Array
31	Invalid Select String Index
32	Requested Select String is empty
33	Select String is too long
34	Invalid Batch Name
35	Object is not on expected Position
36	A200S Protocol non Decimal Digit Received
37	A200S Protocol invalid Termination Character
38	Unknown Object Class
39	No Index found for requested SSL_String
40	Command Value out of Range
41	Parameter Value out of Range
42	HostComm Parser found Invalid Keyword
43	Invalid StrToNum Conversion
44	Parser missing Parameters
45	Invalid Character in Numeric Parameter
46	Parameter Value out of Range
47	Invalid Baud rate Selected
48	PAL Status is < > PAL_IDLE
49	Invalid Sample Number From Remote
50	Answer Count Plunger Move Detection not valid
51	Invalid Plunger Move Detection Parameter
52	Answer Count Needle Magnet not valid
53	Invalid Needle Magnet Parameter
54	Answer Count Head Tool ID not valid
55	Invalid Head Tool Parameter
56	Answer Count Head Board Revision Level invalid
57	Invalid Head Board Revision Level Parameter
58	Answer Count Heater Power Count Invalid

**Table 46.** PAL firmware error codes (Sheet 3 of 6)

<b>Error code number</b>	<b>Error message</b>
59	Invalid Heater Power Parameter
60	Load Firmware not successful
61	RS232 Communication Parameter Error
62	RS232 Channel already Open
63	RS232 Channel not Open
64	RS232 Channel Timeout While Reading/Writing
65	RS232 Parity Error
66	RS232 Framing Error
67	RS232 Overrun Error
68	RS232 RX fifo overflow
69	No serial Channel defined for requested Function
70	Unknown Data Type detected
71	Unknown Item Element detected
72	Unknown Temp Channel
73	Unknown RPM Channel
74	Temperature out of Range
75	Invalid Cycle Detected
76	Answer Count Heater GetRawTemp invalid
77	Invalid Heater GetRawTemp Parameter
78	Parenthesis "(" expected
79	Parenthesis ")" expected
80	Current is too high
81	Speed is too high
82	Acceleration is too high
83	Target position is below low_limit_steps
84	Target position is above high_limit_steps
85	No Syringe Found
86	Syringe Volume out of Range
87	Tray Index out of Range
88	Not Appropriate Object Class
89	Temperature is out of Heater Limits



**Table 46.** PAL firmware error codes (Sheet 4 of 6)

<b>Error code number</b>	<b>Error message</b>
90	Motor Axis Target Position Error
91	Motor Current Fault Phase A
92	Motor Current Fault Phase B
93	Motor Current Fault Both Phases
94	PAL is in PAL_ERROR Status
95	Syringe ID Mismatches
96	Timeout when waiting for answer
97	Answer to big, buffer overflow
98	Timeout during send, SW error
99	Checksum error in received frame
100	Wrong toggle bit in rx frame
101	Invalid procedure parameter >127
102	Could not load head firmware
103	Motion Fault
104	Unknown Atom
105	Not Implemented
106	Object Name is too long, max. 8 Characters
107	Object Named 'None' requested
108	No Syringe installed
109	Invalid SoftTimer
110	SoftTimer Timeout occurred
111	ReportLog Index out of Bounds
112	Invalid Vial Transport mode
113	Invalid Barcode mode
114	Memory Full
115	Found No Vial
116	Address Error detected
117	Bus Error detected
118	Zero Division detected
119	Spurious Interrupt detected
120	Format Error and uninitialized Interrupt

**Table 46.** PAL firmware error codes (Sheet 5 of 6)

<b>Error code number</b>	<b>Error message</b>
121	Input Voltage below trip point
122	Unknown Motor
123	Temperature Sensor failed
124	Valve Position Error
125	Restricted Object Class
126	Heaters are not supported by this Hardware
127	Invalid Host Frame State detected
128	Invalid Host Sequence Number detected
129	Invalid Host Frame Data detected
130	Host Frame ETX Expected
131	Is Not a Hexadecimal Digit
132	Invalid Host Frame Length detected
133	Invalid Host Checksum detected
134	Unexpected STX detected
135	Unexpected ETX detected
136	Unexpected Character detected
137	Needle Guide not Released
138	Wrong Dilutor Syringe inserted
139	Axis stuck on Limit Switch
140	Axis could not detect Limit Switch
141	Axis Low Limit Error
142	Axis High Limit Error
143	Axis Target Position Error
144	Motor Current fault
145	Time Table Full
146	No Tray found at this position
147	z-Axis Collision Error
148	Mandatory Parameter Missing
149	Invalid Vici Valve Drive
150	Invalid Vici Valve Position
151	Invalid Vici Valve Status

**Table 46.** PAL firmware error codes (Sheet 6 of 6)

Error code number	Error message
152	Aux Motors are not supported by this Hardware
153	OutExpBox is currently disabled
154	Output currently used by OutExpBox
155	Invalid Dilutor Motor Status
156	Serial Valves are currently disabled
157	Serial Valve Communication Error
158	Invalid Object Name
159	Invalid Tray Type Group
160	Invalid manual syringe selection
161	Stop on XY Position Error
162	Object used by other object
163	Plunger Position Error Detected
164	Duplicate syringe ID found
165	Washstation is not of type DLW
-1	Unknown Exception received

## External Connectors for Accela Open Autosampler

This section contains information about external connectors for the Accela Open Autosampler.

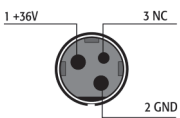
- [Connector Power](#)
- [Connector Fuse](#)
- [Connector LAN](#)
- [Connector AUX1](#)
- [Connector AUX2](#)
- [Connector MODBUS](#)
- [Connector INTERFACE 1](#)
- [Connector INTERFACE 2](#)
- [Connector SER1](#)

## A Reference Information

External Connectors for Accela Open Autosampler

- [Connector SER2](#)
- [Connector TERMINAL or SER3](#)
- [Connector wash station](#)

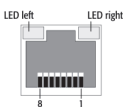
### Connector Power

Pin	Signal name
	
1	+36 V
2	GND
3	NC

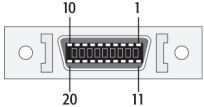
### Connector Fuse

Fuse type	Description
Fuse Type	FST 5 × 20
Rating	6.3 A
Rated Voltage	250 Vac

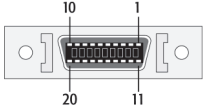
### Connector LAN

Pin	Signal name	DIR	Primary function
			
1	TX+	Out	Transmit Data +
2	TX –	Out	Transmit Data –
3	RX+	In	Receive Data +
4, 5	Not used		Terminated
6	RX –	In	Transmit Data –
7, 8	Not used		Terminated

### Connector AUX1

Pin	Signal name
	
1,2	Motor A1
3,4	Motor B1
6	Temp +
7	Sens
8,9	Heater
11,12	Motor A2
13,14	Motor B2
16	Temp –
17	5 V
15,18,19	GND
20,10	36 V

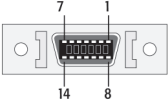
### Connector AUX2

Pin	Signal name
	
1,2	Motor A1
3,4	Motor B1
6	Temp +
7	Sens
8,9	Heater
11,12	Motor A2
13,14	Motor B2
16	Temp –
17	5 V
15,18,19	GND
20,10	36 V

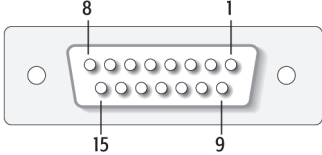
## A Reference Information

External Connectors for Accela Open Autosampler

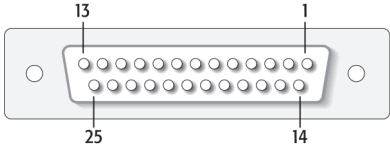
### Connector MODBUS

Pin	Signal name
	
1	TXD
2 – 7	36 V
8	RXD
9 – 14	GND

### Connector INTERFACE 1

Pin	Signal name
	
1	Pwr-Out1 +
2	Pwr-Out2 +
3	SW-Out1 NO
4	SW-Out1 COM
5	SW-Out2 NO
6	Opto-In1 +
7	TTL-In1
8	GND
9	Pwr-Out1 –
10	Pwr-Out2 –
11	TTL-In2
12	SW-Out2 COM
13	TTL-In3
14	Opto-In1 –
15	5 V

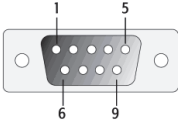
## Connector INTERFACE 2

Pin	Signal name
	
1	TTL-In1
2	TTL-In2
3	TTL-In3
4	TTL-Out1
5	TTL-Out2
6	TTL-Out3
7	Opto-In1 +
8	Opto-In2 +
9	SW-Out1 NO
10	SW-Out2 NO
11	Pwr-Out1 +
12	Pwr-Out2 +
13	5V
14-19	GND
20	Opto-In1 –
21	Opto-In2 –
22	SW-Out1 COM
23	SW-Out2 COM
24	Pwr-Out1 –
25	Pwr-Out2 –

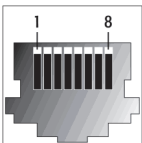
## A Reference Information

External Connectors for Accela Open Autosampler

### Connector SER1

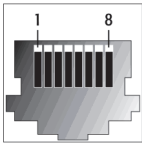
Pin	Signal name
	
1	NC
2	RXD
3	TXD
4	DTR bridged with Pin 7. Special grounding; do not alter.
5	GND
6	NC
7	RTS bridged with Pin 4; do not alter.
8	NC
9	NC

### Connector SER2

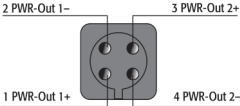
Pin	Signal name
	
1	GND
2	RXD
3	TXD
4	NC
5	NC
6	5 V
7	NC
8	GND



### Connector TERMINAL or SER3

Pin	Signal name
	
1	GND
2	RXD
3	TXD
4	NC
5	NC
6	5 V
7	NC
8	GND

### Connector wash station

Pin	Signal name
	
1	PWR – Out1 +
2	PWR – Out1 –
3	PWR – Out2 +
4	PWR – Out2 –

## System Limitations

This section contains information about system limitations.

### Accela Open Autosampler

You can operate the DLW Option only with an Accela Open Autosampler, which requires PAL Firmware level 4.1.x and the APR Control-xt board.

### Sample Volume

The current version of the DLW Option can handle sample volumes of 2–100  $\mu\text{L}$ .

### Wetted Parts and Material

Some components are not compatible with certain solvents or chemicals that are used in given applications.

### Tubing Internal Diameter (ID) and System Backpressure

Any tubing and component in the flow path, such as an injection valve, contributes to the system backpressure. For example, short tubing with an 0.13 mm (0.005 in.) ID can build up high backpressure when using standard wash pump flow rates.

The wash solvent pumps operate in the range of mL/min. Such a flow rate combined with a high system backpressure caused by small ID tubing will slow down the pump and, as a result, the flow rate.

The described situation is only critical if you select an HPLC setup dedicated for nano flow.

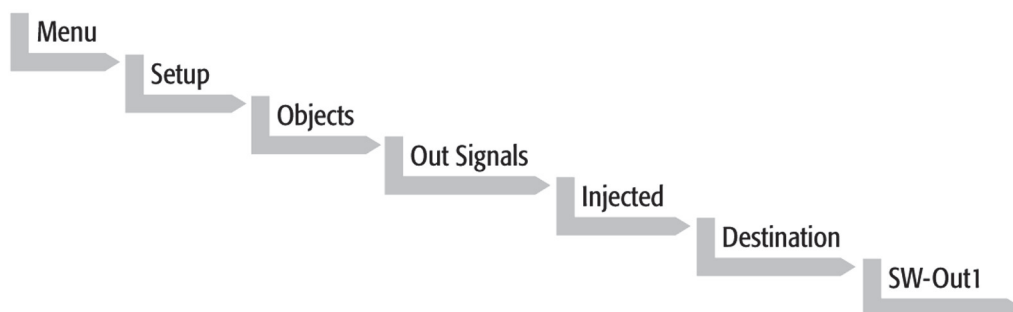
## Interfacing the Open Autosampler to Other Devices

This section contains information about interfacing the Open Autosampler to other devices.

### Synchronization and Output Signals

Synchronization Signals (Sync Signals) are inputs that tell the Accela Open Autosampler when to wait or proceed with a sample-processing step. Output Signals (Out Signals) are sent from the PAL to external devices to indicate status or completion of particular processing steps. These signals are classified as Objects. Physical Events (for example, TTL-In1) are also Objects and may be assigned to named signals. PAL Cycles require that certain signals, such as Start, Inject, and Injected, be defined. For the PAL Events and signal assignments associated with the standard PAL LC-Inj and LC-Cut cycles, see [Table 47](#).

Certain types of integration of the PAL System into data handling or control software, such as ChemStation, Xcalibur, Galaxie or Chromeleon, accomplish synchronization of the Ready/Start signal directly via software control (RS232/LAN).



To assign a physical Event (for example, SW-Out 1) to an existing signal Object (Injected), complete the following menu selections:

The Accela Open Autosampler is shipped with all cycle events predefined, as shown in [Table 47](#). If you need a different physical signal, then you must assign a new Event to the Object signal.

For a signal to be ignored, its corresponding Event must be set to Immediate.

**Table 47.** PAL Events and signal assignments for the standard PAL LC-Inj and LC-Cut cycles (Sheet 1 of 2)

Standard assignment for Open Autosampler injection cycles		Description	Default events	INTERFACE1 pin #
Start JobQueue	(Sync Signal)	Start a Job Queue (Job). Prerequisite is a defined Job with a method assigned. Useful for automated, unattended, timed sampling.	Immediate	
Start	(Sync Signal)	Start a cycle. (Input from GC or data system)	Immediate	
Start2	(Sync Signal)	Continue the cycle GC-Dual for the second sample.	Immediate (Ignore)	
Inject	(Sync Signal)	Inject. READY to INJECT	Immediate (Ignore)	
Inject2	(Sync Signal)	Inject the second sample for GC-Dual cycle.	TTL1 Interface pin7, pin8 (GND)	
Injected	(Out Signal)	Activated at the moment the sample has been injected.	SW-Out1	3 4

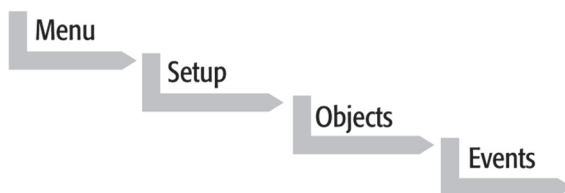
## A Reference Information

Interfacing the Open Autosampler to Other Devices

**Table 47.** PAL Events and signal assignments for the standard PAL LC-Inj and LC-Cut cycles (Sheet 2 of 2)

Standard assignment for Open Autosampler injection cycles	Description	Default events	INTERFACE1 pin #
Injected2 (Out Signal)	Activated at the moment the second sample has been injected.	SW-Out2 Interface pin5, pin12	
Running (Out Signal)	Active as long as the Job Queue is being processed and the PAL is not in an error state.	Off	

If you assign an item of the class Sync Signal to an output signal, such as TTL-In1, then you must define this TTL contact as either Active High or Active Low, as expected by the HPLC system. This is defined in the PAL Firmware class Events.



Select the corresponding item and set the mode accordingly.

If you ordered a dedicated cable, the setting is provided with the schematic description of the LC Sync Cable.

An LC Sync Cable is supplied, which has the PAL Interface 1 connector mounted. The HPLC system side has a 25-pin connector for 600 pump and 1250 pump, and a 2-wire connector for the detectors.

# PAL Loader

This appendix describes how to use the PAL Loader software version 2.5.0 or later.

## Contents

- [Introduction](#)
- [New Features of PAL Loader Software 2.1.x](#)
- [Installation of PAL Loader Software](#)
- [Multiple Shortcuts Assignment](#)
- [PAL Loader Software Operation](#)
- [Troubleshooting](#)
- [Limitations](#)

## Introduction

The purpose of the PAL Loader software is to create a backup file for the complete Accela Open Autosampler firmware. The core software (firmware) and all PAL Firmware Objects, including all settings (Items), local methods, and jobs, are saved as one file.

The PAL Loader software has a second function of loading the firmware or a complete backup file to the Accela Open Autosampler. The PAL Loader application has always been part of the Accela Open Autosampler.

The following PAL Loader software versions are in use:

- PAL Loader version 1.0.1
- PAL Loader version 1.1.1

A new PAL Loader software, version 2.1.0 or later, has been developed for the Accela Open Autosampler.

The Accela Open Autosampler requires PAL Firmware 4.1.x or later, which calls for the new PAL Loader software.

The new PAL Loader software version 2.1.0 (or later) is backwards compatible. The software can be used for all PAL Firmware levels.

The name of the .exe file has been changed intentionally to avoid conflicts when old and new versions are installed on the same computer:

PAL Loader version 1.1.1: PALLOAD.exe  
PAL Firmware level up to and including 3.x.x

PAL Loader version 2.1.x: Setup PALloader.exe  
PAL Firmware level as of 4.1.x or later  
but backwards compatible with earlier versions.

## New Features of PAL Loader Software 2.1.x

The following features, new since version 1.1.1, are available in the new PAL Loader software:

- The installation software is executable directly from a CD-ROM or any transportable storage device, such as a memory stick.
- The software installation procedure is different from that of Loader version 1.1.1. If they do not already exist, directories are created in the Microsoft User Profile.
- The read/write speed for a file is significantly higher. The initial COM Port settings, such as the baud rate, are identical to previous versions. The baud rate is adjusted dynamically during program startup and is displayed on the status line, in the lower right corner of the application window.
- You can create multiple shortcuts and for each one assign a dedicated IP or COM address. This setup facilitates the handling of a number of PAL Systems from a single computer.

## Installation of PAL Loader Software

This section contains information about installing the PAL Loader software.

### Requirements

PAL Loader software version 2.1.x requires Windows 2000, XP, or Vista™. It does not support earlier Windows versions, such as Windows 95, 98, or NT. If installation is necessary in this environment, use PAL Loader version 1.1.1.

Software communication from a computer to the PAL System can be done one of two ways:

(Serial control, requiring a serial cable provided with the PAL System, crossed wiring)

- P/N 00950-01-00332 (cable length 3 m, standard cable)

–or–

- PN: 00950-01-00333 (cable length 6 m, optional cable length)

## PAL Loader Software Version 2.1.x Installation

The software is provided with every PAL-xt System, with an enclosed CD-ROM.

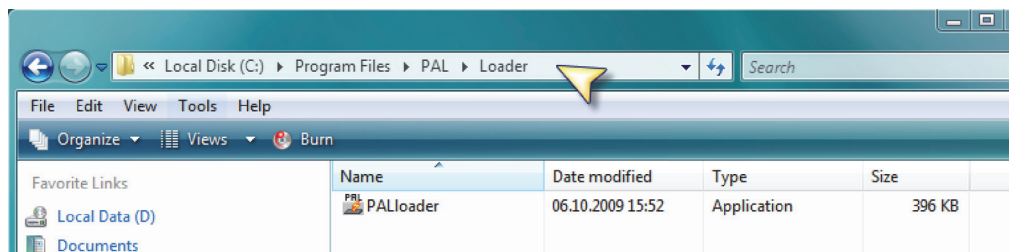
The PAL Loader.exe file is embedded in a wizard software package. Starting the wizard executes the installation software. If no changes are made to the proposed installation path, the Loader version 2.1.x uses the following default installation path (Figure 74) to install the software:

*drive:\Program Files\PAL Loader*

**Note** You can start the wizard for the installation software by double-clicking the icon of the file Setup PALloader.exe.

You can use the second file, PALloader.msi, for the installation as well. Double-click the icon of the file, right-click, and choose Install from the shortcut menu. This path for installation requires an up-to-date Windows Installer. If the current version is not installed, a corresponding error message appears.

**Figure 74.** Installation path for the PAL Loader software

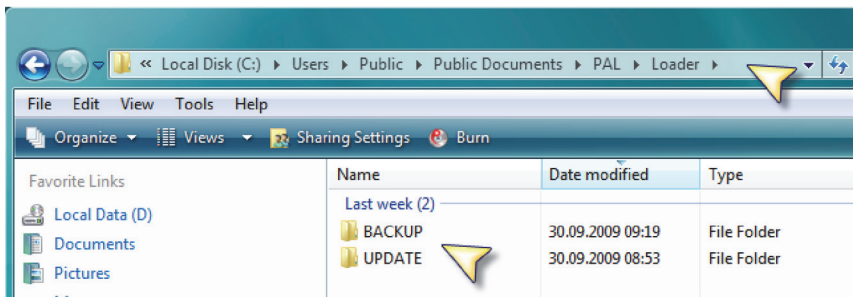


The two subfolders Backup and Update are automatically created at installation in this directory:

*drive:\Users\Public\Public Documents\PAL\Loader*

The directory allows writing temporary files, with no write protection and, in a network environment, the User Profile is always provided to the user at logon.

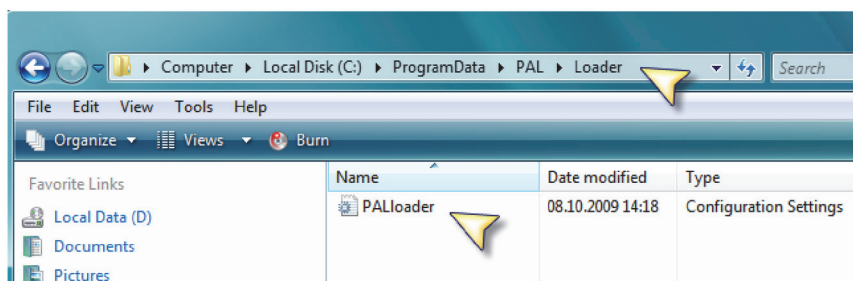
**Figure 75.** Installation path for public documents



A Loader INI file is created at installation and stored in this directory:

*drive:\ProgramData\PAL\Loader*

**Figure 76.** Installation path for PAL Loader INI file



Note that the folder ProgramData can be hidden if a standard installation of the Windows Operating system is executed.

## Multiple Shortcuts Assignment

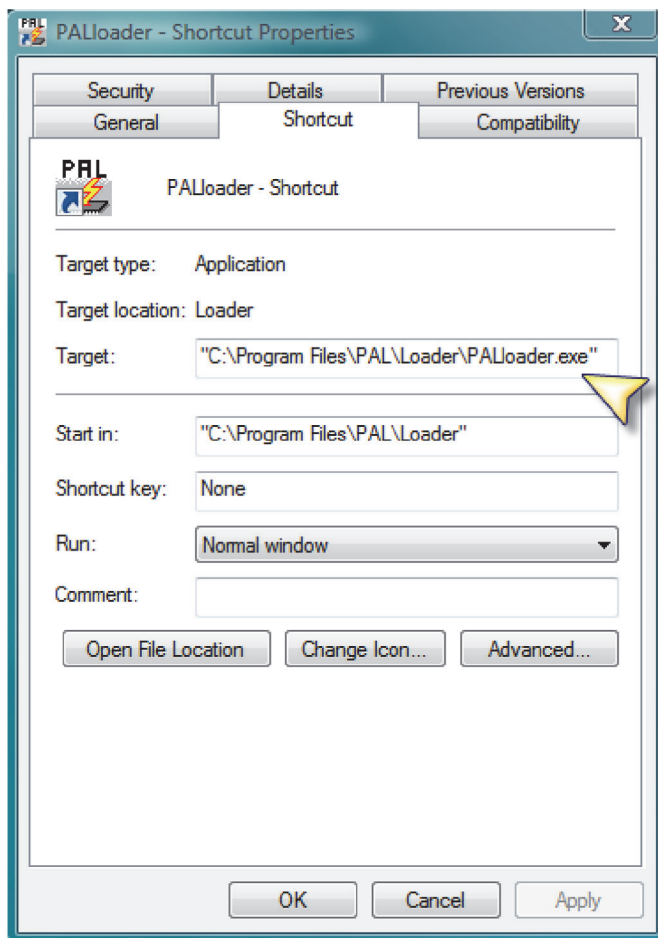
The PAL Loader version 2.1.x or later provides for the handling of multiple shortcuts on the desktop; each one can be assigned to a specific TCP/IP address or COM port. A mix of the two communication protocols is possible as well.

This setup facilitates the handling of a number of PAL Systems from a single computer.

You can create a shortcut of the PALloader.exe file to the desktop. After you create the shortcut, right-click the shortcut icon and choose Properties from the shortcut menu. See [Figure 77](#).



**Figure 77.** Shortcut Properties dialog box



The Target command line uses the installation path of the PAL Loader software:

*drive:\Program Files\PAL\Loader\PALloader.exe*

**Tip** Create a shortcut from the data application file located in the above described path (in Explorer) and not from the PAL Loader desktop icon.

## Dedicated Shortcut for Serial Communication

Add this extension to the command line:

`... .exe /CommunicationType=0 /ComPort=x`

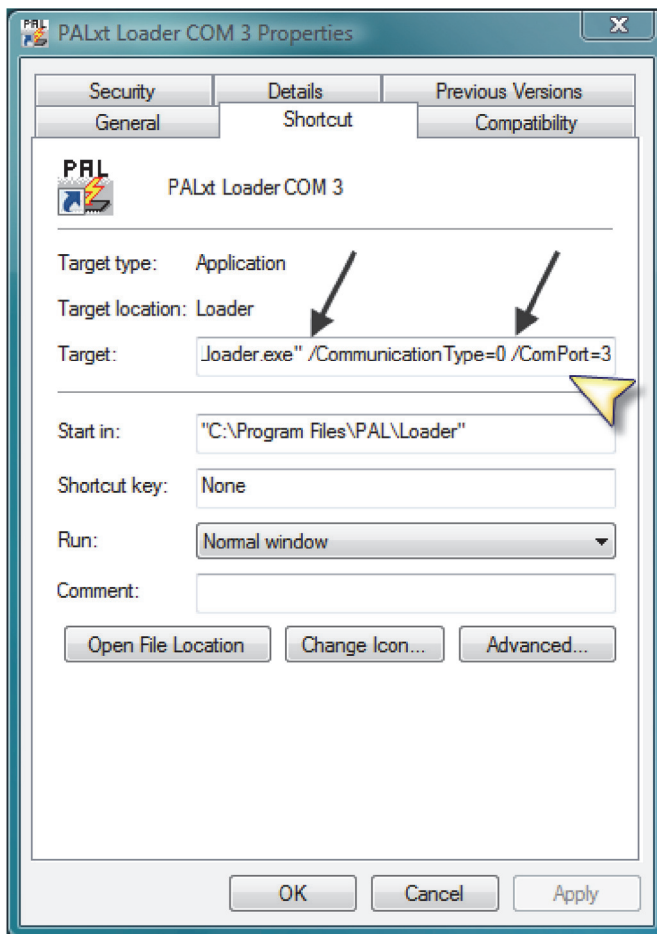
See [Figure 78](#).

Where:

- Type=0: Serial Communication Protocol
- x= Com Port Number

**Tip** After ...exe", insert a blank space and a slash ("/", not a backslash), then ...**Type=0** and another blank space and slash, "/".

**Figure 78.** Extending target command for serial communication



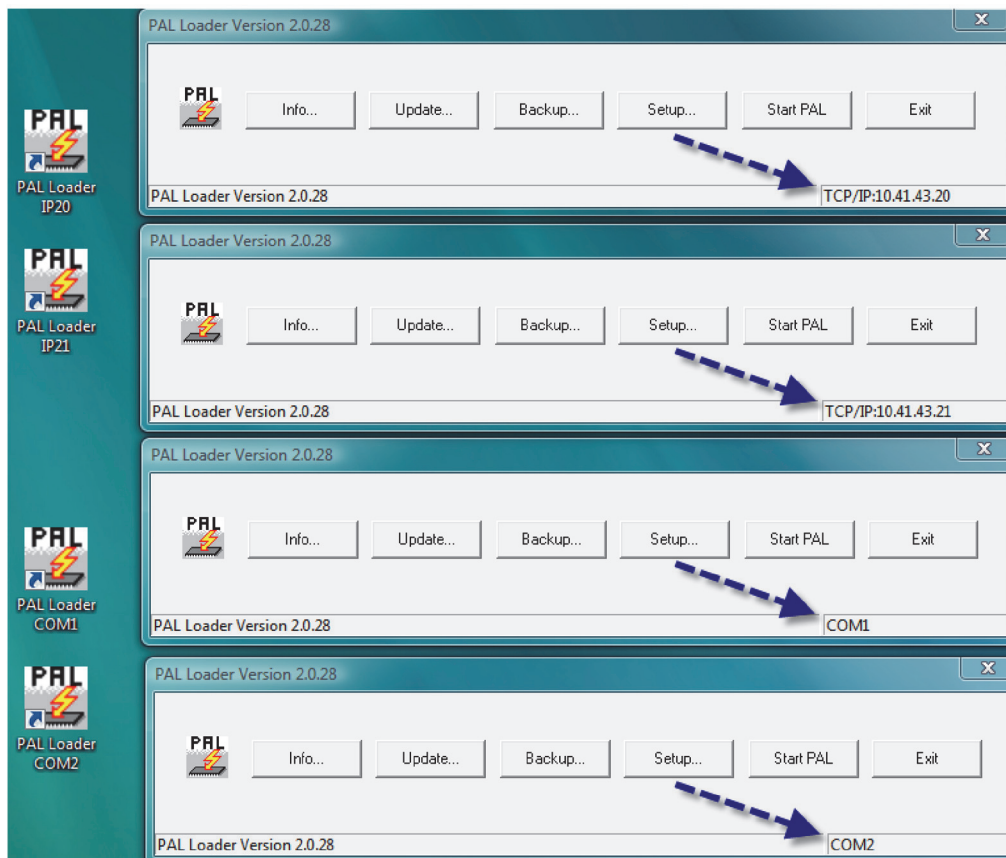
## Creating Multiple Shortcuts

You can use the same procedure as described in the previous section to create and save several shortcuts on the desktop.

For each shortcut icon, right-click and choose Rename from the shortcut menu to define a unique name. [Figure 79](#) shows an example of two LAN shortcut icons with IP address 20 and 21 as well as two serial shortcut icons with COM port numbers 1 and 3.

Double-clicking the icon opens the PAL Loader software with the assigned communication protocol and address (port).

**Figure 79.** Example of multiple shortcuts



If several PAL Loader applications are open, as shown in [Figure 79](#), and you close them successively, the address of the last closed application is written back to the Loader INI file.

Reopening any of the PAL Loader applications with a dedicated communication protocol and address (port) will use the specified setting and overwrite the INI-file settings.

## PAL Loader Software Operation

This section contains information about how to operate the PAL Loader software.

### Starting the PAL Loader Software

Click the PAL Loader icon located on the desktop.

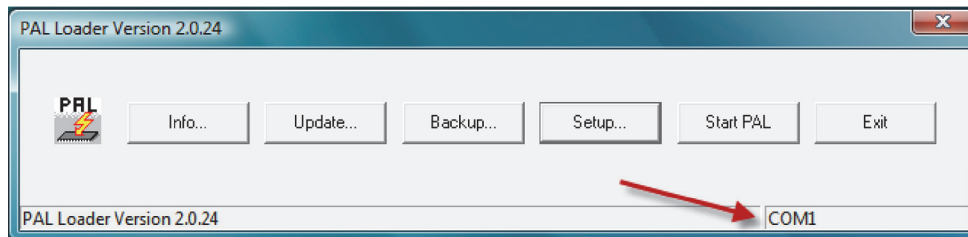
**Figure 80.** PAL Loader software icon



The application opens and displays the various user buttons (Figure 81).

The previously used communication mode is displayed in the status bar.

**Figure 81.** PAL Loader application buttons



## Setting up COM Ports

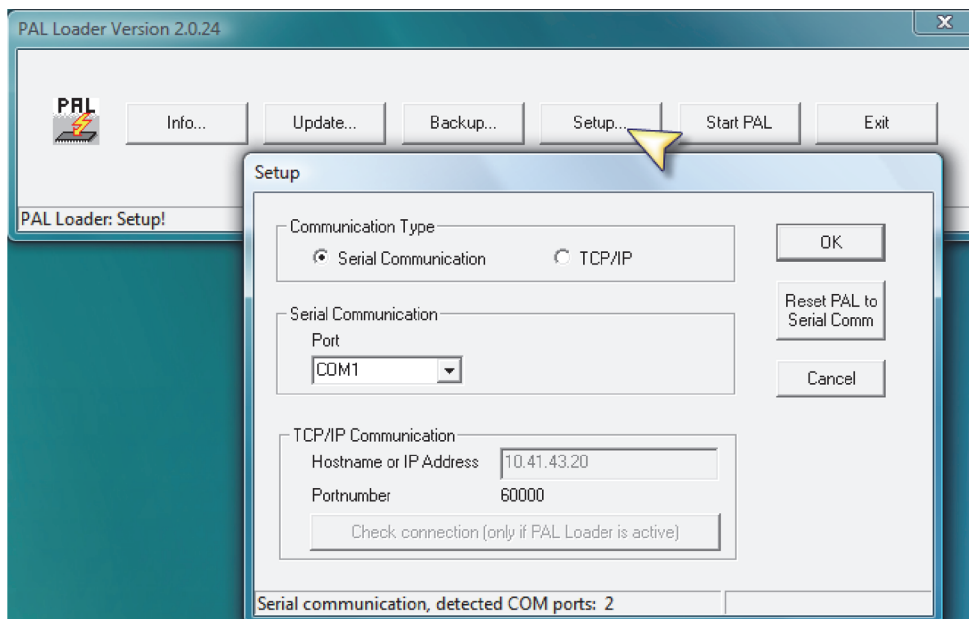
Clicking the Setup button opens the dialog box to select serial communication (Figure 82).

The COM Port is user selectable. The message displays the number of COM ports detected.

**Note** The limitation of PAL Loader software version 1.1.1—that only ports COM1 to COM4 can be addressed—has been eliminated in version 2.1.x.

Also, COM ports do not need to be in consecutive order.

**Figure 82.** Setup dialog box for Serial Communication



## COM Port Settings

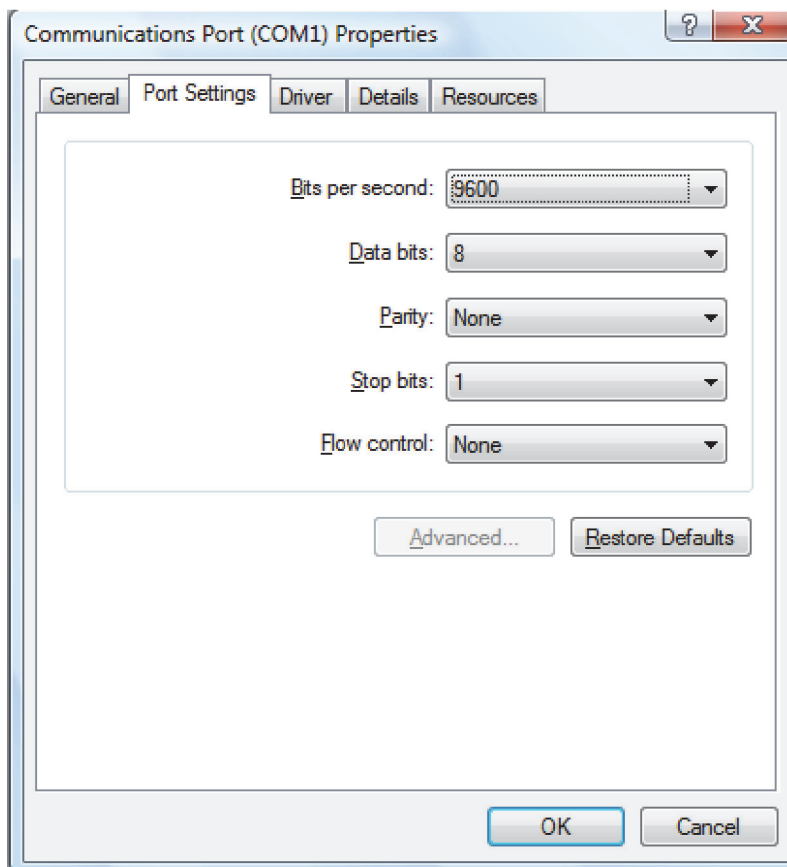
The port settings are normally defined in the following:

Windows\Control Panel\Device Manager\Ports

The standard settings for serial communication are shown in [Figure 83](#).

**Note** The application (PAL Loader software) actively sets the communication parameters, and the baud rate is set for optimized use.

**Figure 83.** COM1 Properties dialog box – Port Settings page



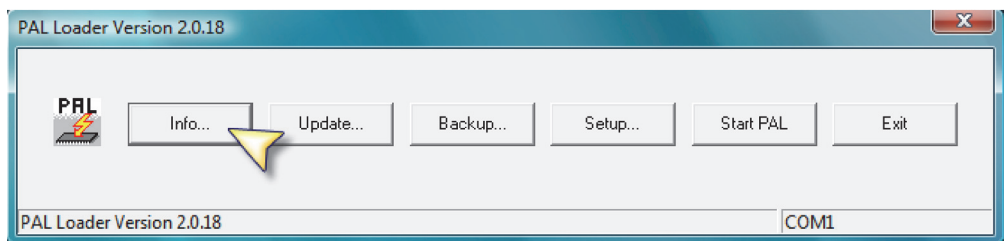
## Serial Communication

This section contains information about serial communication.

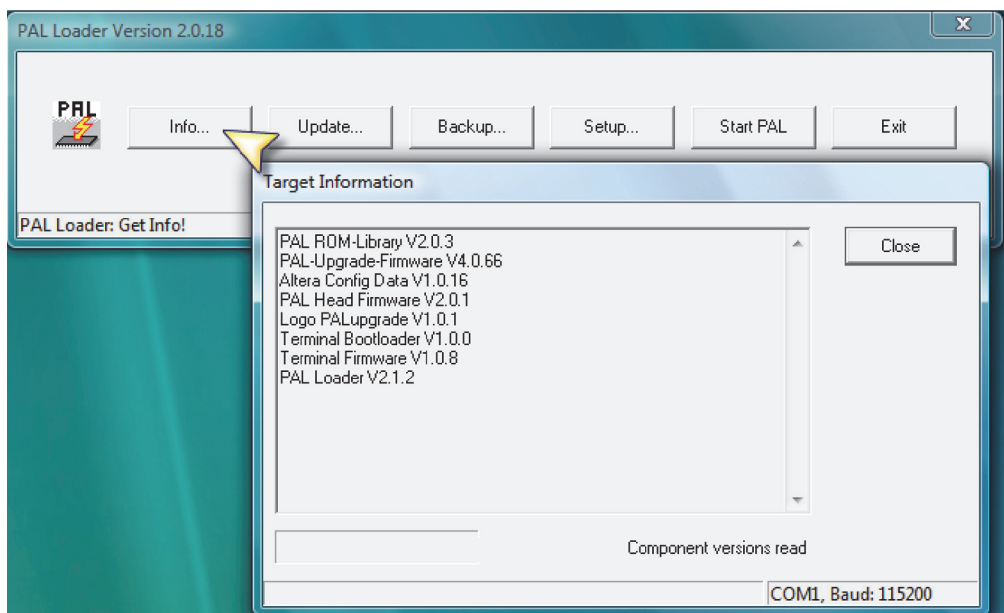
### Establishing Serial Communication

To quickly check whether serial communication with the PAL System can be established, click **Info** (Figure 84 and Figure 85).

**Figure 84.** Clicking Info



**Figure 85.** PAL information received



When the PAL Loader application receives information from the Accela Open Autosampler, the red LED on the back of the electronic board blinks. This blinking LED indicates that the processor is in so-called Loader Mode, which means that the processor enables receiving the firmware or writing to memory. For more details, see [LED Status and Position](#).

**Note** The Target Information list displays the version levels of the different components of the Accela Open Autosampler.

The various types of information received might be important for troubleshooting. For example, if the line for the PAL Firmware release is not shown, the firmware is most likely corrupt.

## LED Status and Position

The following are the various states for the LED indicator on the Open Autosampler:

- Turned off: The autosampler is in normal operating mode.
- Blinking or flashing: The autosampler is prepared in Loader Mode to create a backup or update—that is, to receive data from, or write to, memory.
- Constantly on: Data transfer process is in progress.

The position of the LED in the APR Control-xt board is between the buzzer and the connector for the wash station.

## Explanation for the Communication Protocol Settings

**Table 48.** Accela Open Autosampler communication parameters

Parameter	Explanation	Remarks
CommPort	Setup for serial (SER1) or LAN communication	If serial communication is selected, all other parameters in PAL-xt Firmware, Communication class, are ignored.

## Reset PAL-xt to Serial Communication

If the PAL-xt communication is set to LAN and for any reason the PAL Firmware became corrupt, you can reset or force back the communication protocol to Serial by using the button, Reset PAL to Serial Comm, in the Setup dialog box of the PAL Loader ([Figure 82](#) on [page 162](#)).

## Backing Up the PAL Firmware

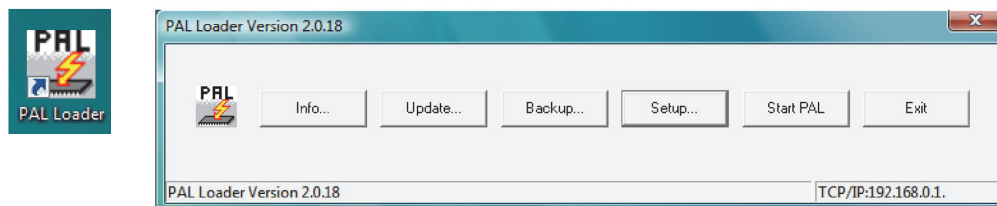
**Note** Perform the following backup procedure after completing installation of a new PAL System. In addition, periodic system backups ensure that a PAL updated configuration file (PAL Firmware backup) is available at any time.

**IMPORTANT** Do not move any part of the PAL System, such as the *y-z*-axis assembly and, if possible, do not use other computer functionality during the backup and update process. This will cause delays and might freeze the window.

### ❖ To create a backup file

1. Shut off power to the PAL System.
2. Connect the cable corresponding to the communication mode.
  - Serial cable (RS232) between PAL SER1 and the computer Serial port (COM)  
PN: 00950-01-00332 (standard cable provided with PAL System) or  
PN: 00950-01-00333 (optional, cable length 6 m)
3. Switch on the PAL System power.
4. Double-click the **PAL Loader** icon (Figure 86). The PAL Loader startup window appears.

**Figure 86.** PAL Loader icon and startup window



5. Click **Backup**. The Backup Target Memory dialog box opens (Figure 87).

The path (storage location) and the file name to be assigned to the new backup file appear.

You may customize the path, file name, or both.

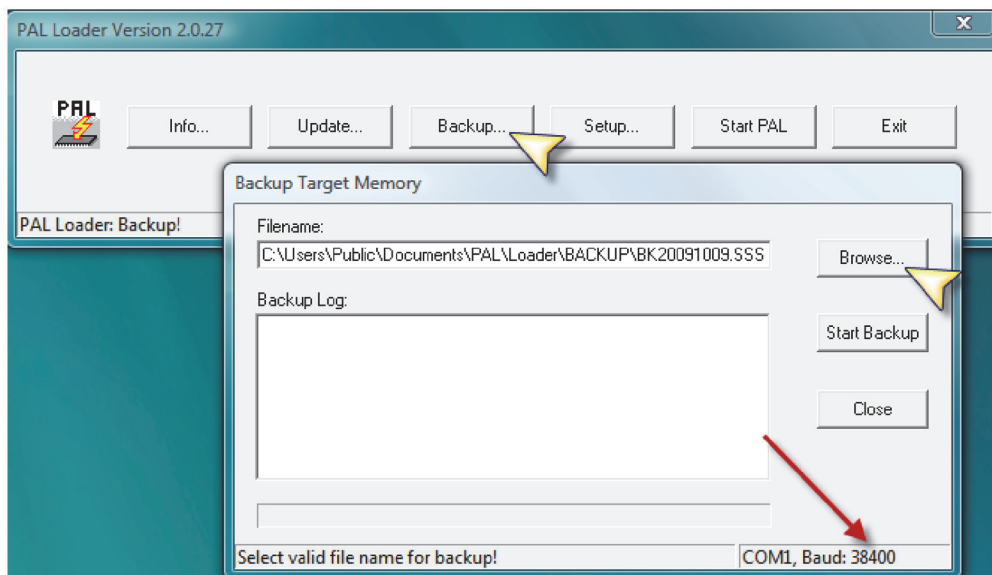
Default file name: BKdate.sss (date: *YYMMDD*)

**Note** You can also use the Info button to establish and check communication with the PAL System using the PAL Loader software.

If you cannot establish communication with the PAL System, see “[Troubleshooting](#)” on page 85.



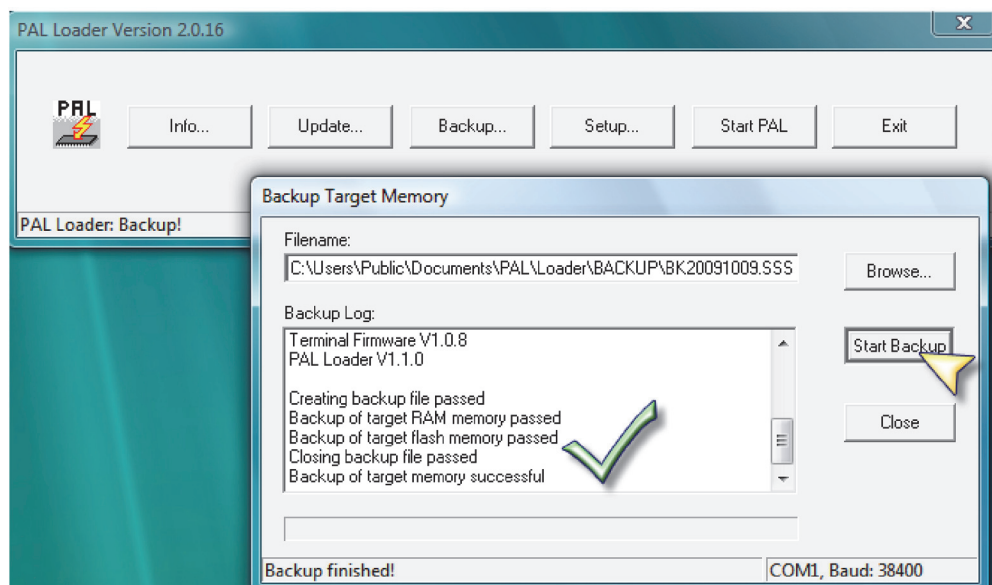
**Figure 87.** Starting the backup process with serial communication type



**Note** The displayed baud rate in this example of 38400 is the baud rate last used. Do not confuse this with the COM port setting, which is a baud rate of 9600. The application optimizes the speed, depending on work load and processor capabilities. The dynamic increase in the baud rate can reach as high as 115 200. The last used baud rate writes back to the PAL Loader INI file. Opening this window from the PAL Loader application displays this last used rate.

6. Activate the process by clicking **Start Backup** (Figure 88). The process is initialized and takes approximately 3–5 minutes.

**Figure 88.** Backup process for serial communication completed



A backup file is created, the RAM and flash memory are backed up, and the newly created file is closed. After the process is complete, the message: Backup of target memory successful is displayed.

7. If you receive the message, “Backup of target memory not successful,” repeat the last step and check the communication by clicking Info.

## Updating the PAL Firmware

**IMPORTANT** Do not move any part of the Accela Open Autosampler, such as the  $x$  axis, and if possible do not use other computer functions during the backup and update process. This will delay the process and might freeze the application.

Use the following procedure to upgrade the PAL Firmware to another level. Keep in mind that the existing firmware data, such as parameter settings or methods, is overwritten or lost. This is especially important for users who work directly from their local terminal and create methods and jobs within the PAL System. In such cases, ensure that the methods are saved before starting an update procedure.

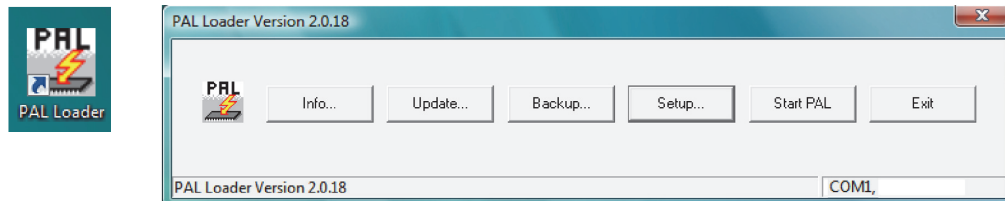
**Note** Best practice dictates creating a backup file as described above before performing an update of the PAL Firmware. This backup file is for backup purposes only. If an important point is overlooked, recovery is possible by reloading the backup file onto the Accela Open Autosampler.

### ❖ To create an update file

1. Shut off power to the Accela Open Autosampler.
2. Connect the cable corresponding to the communication mode:
  - Serial Cable (RS232) between PAL SER1 and the computer Serial port (COM)  
PN: 00950-01-00332 (standard cable provided with PAL System) or  
PN: 00950-01-00333 (optional, cable length 6 m)
3. Double-click the **PAL Loader** icon (Figure 89).

The PAL Loader startup window appears. The previously used communication mode is displayed in the status box.

**Figure 89.** PAL Loader icon and startup window



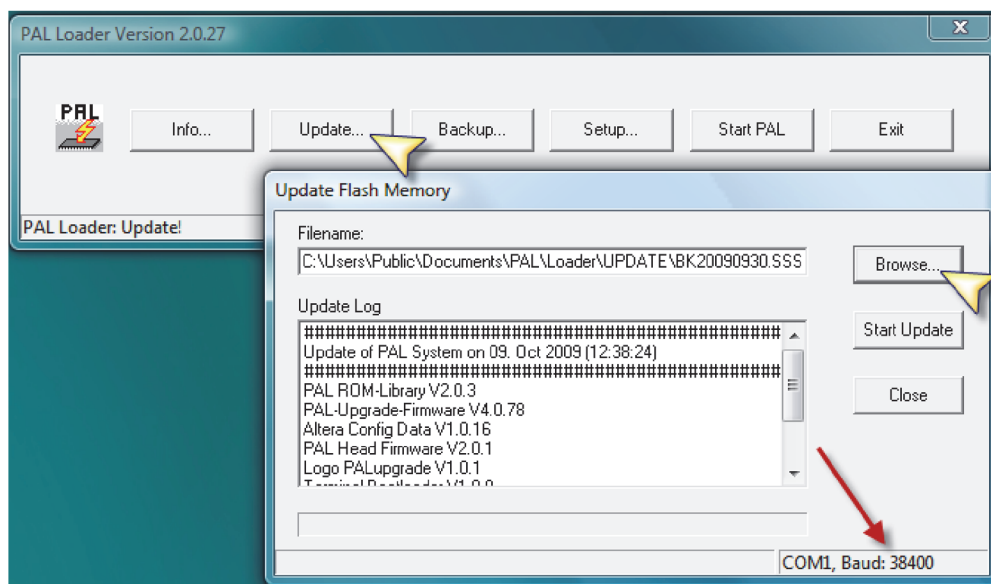
4. Click **Update**.

The Update Flash Memory dialog box opens (Figure 90).

5. Browse to the appropriate file (extension .sss) to upgrade the PAL Firmware.

**Note** To establish and check communication from the Accela Open Autosampler using the PAL Loader software, you can also use the Info button. If communication with the Accela Open Autosampler cannot be established, see “[Troubleshooting](#)” on page 85.

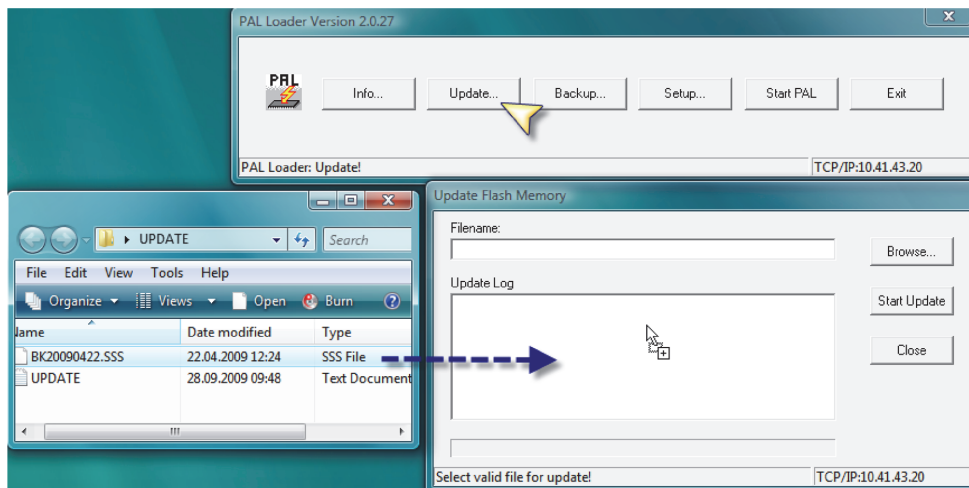
**Figure 90.** Starting the update process with a serial communication type



**Note** The displayed baud rate in this example of 38 400 is the last used baud rate. Do not confuse this with the COM port setting, which is a baud rate of 9600. The application optimizes the speed, depending on work load and capabilities of the processor. The dynamic increase of the baud rate can reach as high as 115 200. The last used baud rate writes back to the PAL Loader INI file. Opening the PAL Loader window displays this last used rate.

6. In PAL Loader software version 2.1.x, load or activate a file (with extension .sss) by dragging it to the window (Figure 91).

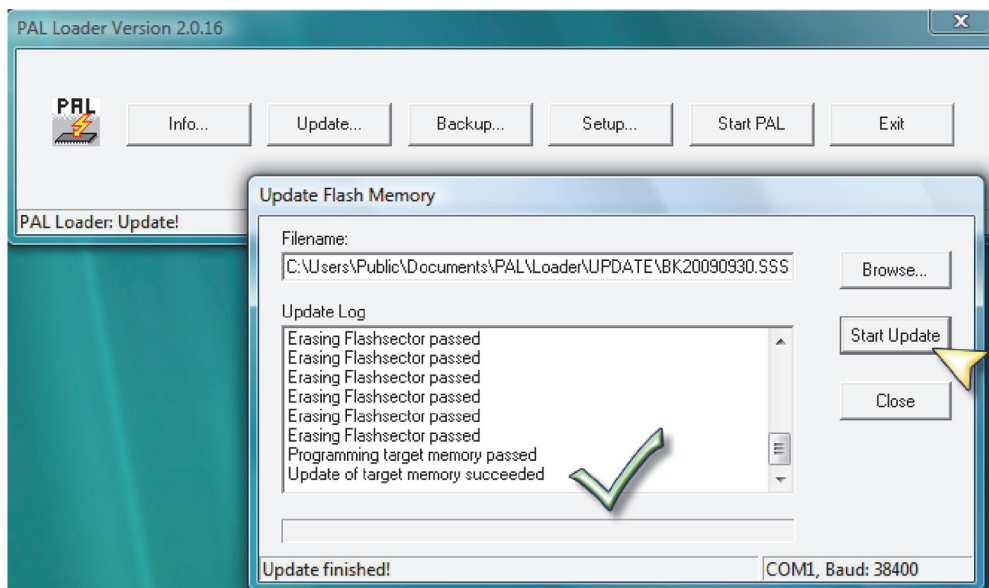
**Figure 91.** Drop a file into the Update window



7. Activate the process by clicking **Start Update** (Figure 92).

The process is initialized and takes approximately 3–5 minutes.

**Figure 92.** Update process for serial communication completed



The RAM and flash memories are erased and replaced by the new data. When the process is complete, this message appears: Update finished (from the task bar).

If there is a problem, a message similar to the following appears: Programming target memory failed at 0x000A2D00.

The Update log entry memory failed #####.

You can try repeating the last step, using the following options:

- Retry with old file.
- Load new file.

You might check the communication by using the Info button.

## Troubleshooting

This section contains information about troubleshooting the PAL Loader software.

### Serial Communication with PAL Loader Software

**Note** You can use PAL Loader software version 2.1.x for the PAL System. It is backwards compatible with PAL Systems.

If you cannot establish serial communication with the PAL Loader software version 2.1.x, check the following points:

- Is the COM port on the computer available and correctly addressed?

Are default settings verified?

- Is the PAL Loader software COM port setting correct?

If no RS232 (serial) port is available on the computer, use a USB/Serial converter. Use the type recommended by CTC Analytics: PNo.: USB-Ser2.

The PAL RS232 cable has a crossed wiring connection.

PN: 00950-01-00332 (3 m length, standard cable) or

PN: 00950-01-00333 (6 m length, optional cable)

### How to Force Loader Mode Manually

If basic conditions are met but you still cannot establish communication from the computer to the Accela Open Autosampler, you must force the PAL System into Loader mode.

This action might be necessary if the PAL Firmware has been partially corrupted.

#### ❖ To force the Loader Mode manually

1. Turn off the power from the PAL System.
2. Turn on the power again.
3. Turn off the power from the PAL System within a time window of approximately 3 to 5 seconds again (turn it off before the startup logo is displayed on the screen).
4. Turn on the power again.

Successful activation of the Loader Mode is indicated by the blinking red LED at the back of the PAL *x* axis.

There are two blinking patterns for the LAN and COM (serial) modes for the Control-xt board only. The LAN mode exhibits a fast, double-blinking signal, whereas the COM (serial) mode shows a pattern with one blinking signal at a constant time interval. See a visual representation of the patterns in [Figure 93](#).

**Figure 93.** LED flashing pattern for LAN and COM modes



Double blinking - LAN Mode



Single blinking - COM (Serial) Mode

Location of LED:

- For PAL Systems, APR CPU or APR Control boards:  
LED is beside the DC connector.
- For the PAL-xt System, APR Control-xt board:  
LED is between the buzzer and wash station.

## Reset PAL to Serial Communication

If the PAL-xt communication protocol is set to the LAN type, it is possible that the link between the PAL-xt software and the PAL-xt System can no longer be established. One cause for such a failure is a corrupt PAL Firmware.

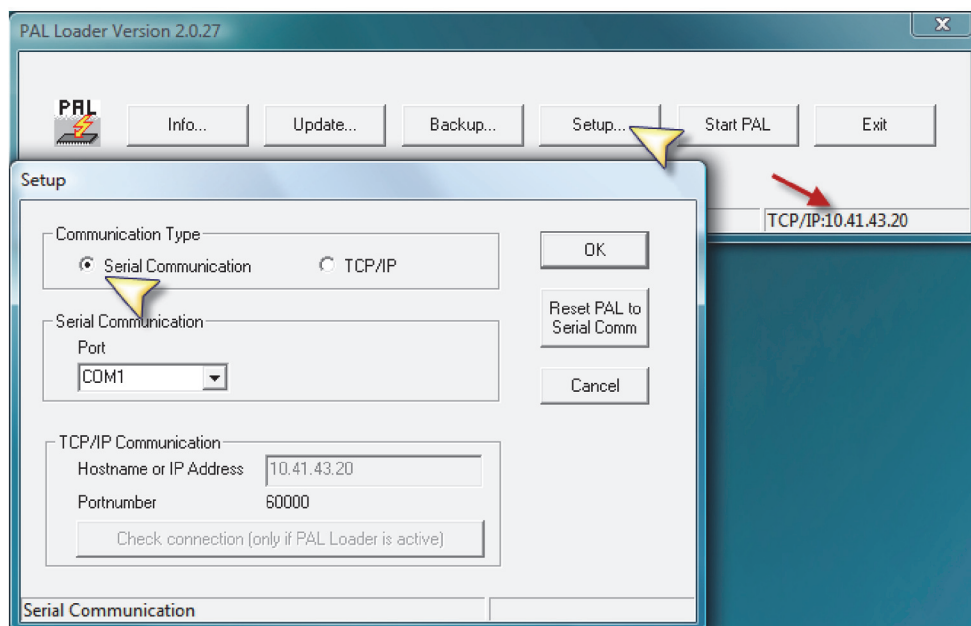
Under these circumstances, you must be able to execute an “emergency” command. This function is set in the PAL Loader software Setup dialog box using the Reset PAL to Serial Comm button.

## How to Reset the PAL to Serial Communication

### ❖ To reset the PAL to serial communication

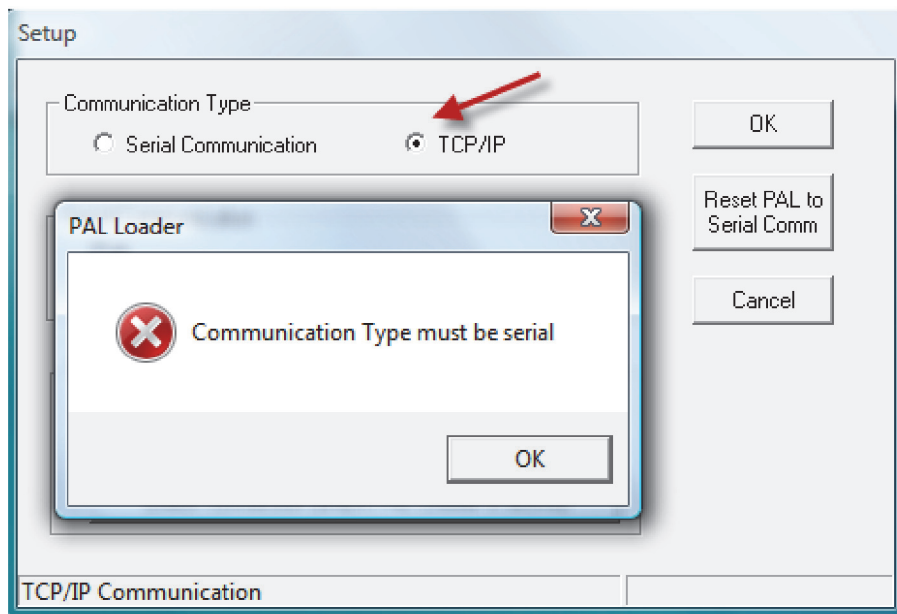
1. Start the PAL Loader application (version 2.1.x or later) by double-clicking the corresponding icon.
2. Click **Setup** and change the Communication Type from the TCP/IP to the **Serial Communication** option (Figure 94).

**Figure 94.** Activating Setup for resetting the PAL-xt



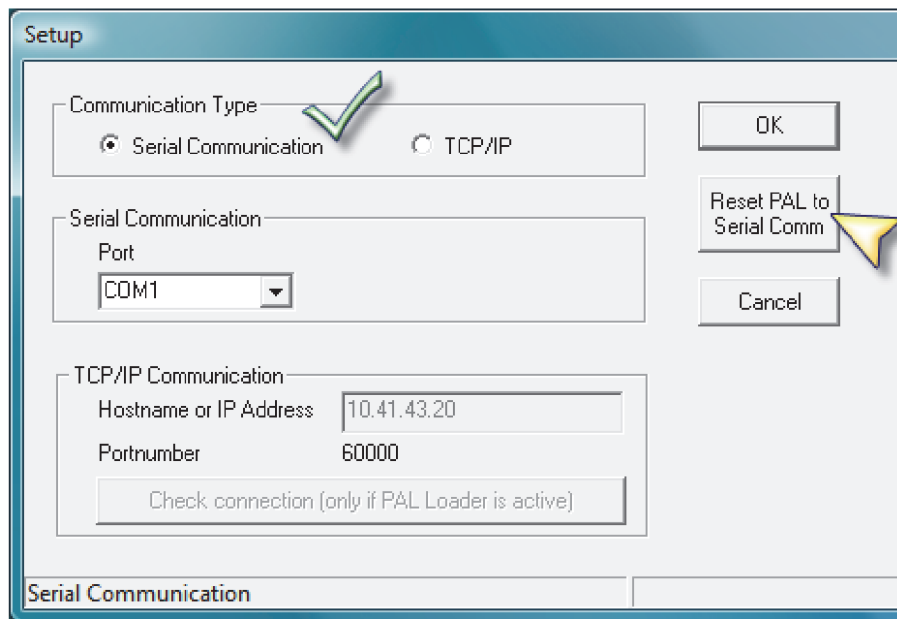
If the Communication Type is not set to Serial, an error message appears (Figure 95).

**Figure 95.** Preparing to reset the PAL-xt



3. After selecting the Serial Communication Type option, click **Reset PAL to Serial Comm.**

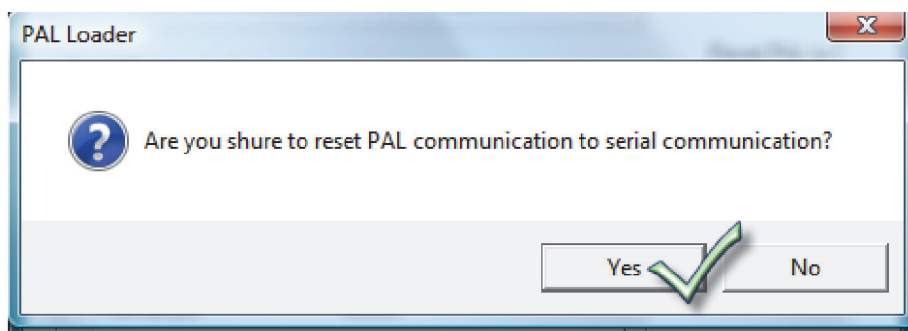
**Figure 96.** Activating Reset PAL to Serial Communication





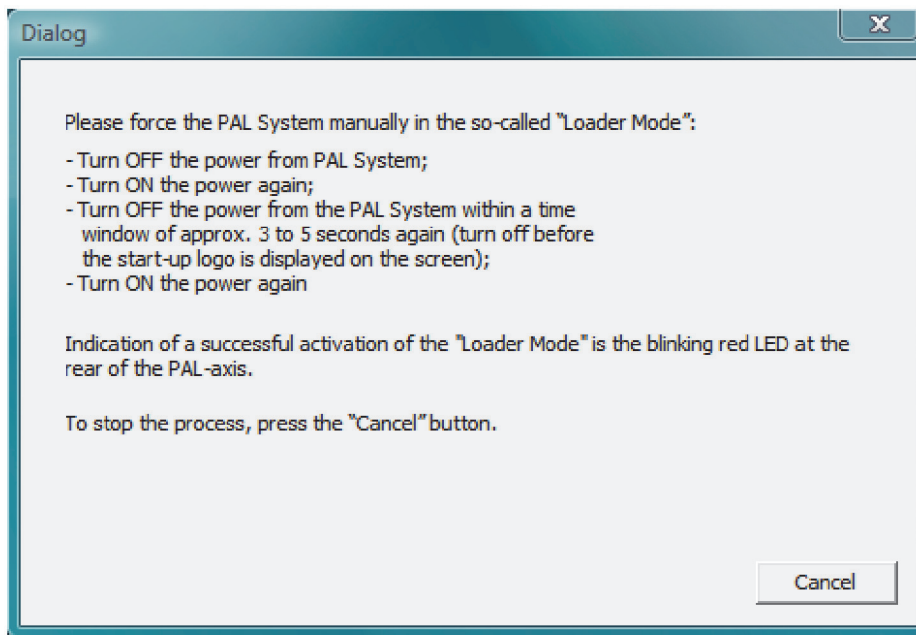
A confirmation box opens (Figure 97).

**Figure 97.** Confirmation for the reset process



4. Click **Yes** to continue the reset process. A second window appears with instructions to manually turn off and on the power of the PAL-xt System (Figure 98).

**Figure 98.** Instructions to force the PAL-xt System into Loader Mode



5. Click **Cancel**.

After successful activation of the Loader Mode, the communication type changes automatically to Serial Communication. A beep tone signals the completion of the reset procedure.

The PAL-xt System is now prepared to perform a PAL Firmware update.

## Erratic Behavior of Software or Flash Memory

In some rare cases the red LED might continue to blink after a PAL Firmware update. You cannot start the PAL System from the PAL Loader Start PAL button or by powering it off/on. If such a case occurs, take the following steps:

1. Power down and wait at least 30 seconds before the unit is powered up again. If the PAL System is not restored, go to the next step.
2. Reload a complete PAL Firmware backup file. A complete backup file means core software including all PAL Firmware Objects data, content of RAM, and flash memory.  
If this action is still unsuccessful, go to the next step.
3. Power down the unit again and remove the board (APR CPU, APR Control, or APR Control -xt). Before you handle an individual board, be careful of static charge.
4. Connect the DC cable, the spiral cable for the terminal, and either the Serial or Ethernet cable to the board on-the-fly.
5. Start the software UPDATE process again.
6. If none of these actions is successful, change the board.

Any severe error requiring changing the board typically means that the Loader section of the flash memory can no longer be addressed, which results in communication failure.

**Note** A PCB, APR CPU, APR Control, or APR Control-xt, which has not been connected to current for a period of six months or longer, can result in a low battery level. Because a low battery will no longer support the RAM memory, data loss is likely.

7. Charge the battery overnight (36 V), and reload a complete PAL Firmware backup file (core software and PAL Firmware Objects).

## Failure to Complete BACKUP or UPDATE Process

Avoid touching the PAL System during the read or write processes (BACKUP or UPDATE). Do not move, for example, the  $y$ -,  $z$ -axis assembly during a process. This will delay its completion and could result in an error message and stopping the process completely.

Furthermore, the window can freeze (shown in a frozen progress bar). This might also happen if the computer processor is very busy with other tasks at the same time. Try to avoid using any other computer function during the process.

## Limitations

Known limitations to consider for troubleshooting purposes relate to the Windows operating system. The PAL Loader software version 2.1.x is compatible with the following systems only:

- XP Professional (Service Pack 2)
- Vista Enterprise (Service Pack 1)
- Windows 7 beta build 7000, preliminary tests fulfilled



## PAL Object Manager

This appendix describes how to install, set up, and use the PAL Object Manager software. This appendix is intended for frequent Accela Open Autosampler users or new users who are experienced at configuring and using automated systems to run existing analytical methods. Before using the PAL Object Manager, make sure you have properly installed and set up the Accela Open Autosampler using the instructions in “[Installing the Accela Open Autosampler](#)” on [page 11](#).

### Contents

- [Installing PAL Object Manager](#)
- [Getting Started with PAL Object Manager](#)

**IMPORTANT** Before using the PAL Object Manager software to make changes to your system, back up the current system definition using the PAL Loader software. If problems occur while using the PAL Object Manager software, repair the system using PAL Loader to restore the back up configuration file. See “[PAL Loader](#)” on [page 155](#).

Because this software can change your system to the point that the autosampler and Thermo PAL does not function properly, make changes to objects only if you are an advanced user and understand how the changes you make can impair system functioning.

Thermo Fisher has defined the hardware and software configuration for your autosampler. Changes to this configuration can impact system functioning.

When you add a new hardware module to the Accela Open Autosampler, use the PAL Object Manager software to load new PAL firmware objects onto the system to support those modules. The Object Manager mode also helps create and maintain object lists from an existing Accela Open Autosampler.

After adding an object to the Accela Open Autosampler, choose **Restart Pal** from the File menu to start the system so other software applications can use the object.

The latest PAL Object Manager software is backwards compatible. You can use the software with all PAL firmware levels.

## Installing PAL Object Manager

Installing PAL Object Manager software requires three steps.

- [Verifying System Requirements](#)
- [Installing the Object Manager Software to the Local Hard Drive](#)
- [Installing the PAL Object Lists](#)

### Verifying System Requirements

Before installing the PAL Object Manager software, your system must meet these minimum requirements (Table 49).

**Table 49.** System requirements for PAL Object Manager

System	Minimum requirements
Software	<p>PAL Object Manager software is compatible with these Windows operating systems:</p> <ul style="list-style-type: none"><li>• Microsoft Windows 2000</li><li>• Microsoft XP Professional (Service Pack 2)</li><li>• Microsoft Vista Enterprise (Service Pack 1)</li><li>• Windows 7 beta build 7000, preliminary tests fulfilled.</li></ul>
Hardware	<p>For serial control communication from the computer to Accela Open Autosampler, use one of these crossed-wiring serial cables provided with the Accela Open Autosampler:</p> <ul style="list-style-type: none"><li>• P/N 00950-01-00332 (cable length, 3 meters)</li><li>• P/N 00950-01-00333 (cable length, 6 meters)</li></ul> <p>If the desktop or laptop computer has no serial port (COM port) installed, use a conversion adapter USB/Serial.</p>

### Installing the Object Manager Software to the Local Hard Drive

The PAL firmware objects are grouped in various folders that must be linked to the Object Manager software. To avoid any conflicts with the link, load the software to your local hard drive before copying the object folders.

You must have administrative rights to install the software.

❖ **To install the Object Manager software to the local hard drive**

1. Put the Object Manager CD in your computer.
2. To start the installation, double-click the **PAL Object Manager setup.exe** file.

The installation wizard begins installing the software.

3. If there is no change to the proposed installation path, install the software to the default location:

C:\Program Files\PAL\Object Manager

## Installing the PAL Object Lists

After installing the Object Manager software, install the PAL Object Lists folder. This folder contains all PAL firmware object lists in alphabetical order by class, for example, syringes, trays, and wash stations.

Install the PAL object lists folder in the same directory as the PAL Object Manager software.

❖ **To install the PAL Object Lists folder**

1. Find the Pal Object Lists folder on the CD.
2. Copy the folder and paste it into the same folder on your hard drive where the PAL Object Manager resides. The default location is C:\Program Files\PAL\Object Manager.

## Getting Started with PAL Object Manager

This section describes how to get started with PAL Object Manager software.

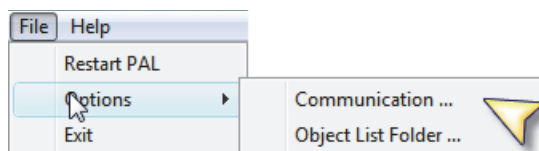
- [Setting Up the Communication Mode](#)
- [Troubleshooting the Serial Communication Mode](#)

## Setting Up the Communication Mode

Use the File menu in the PAL Object Manager window to set the communication mode.

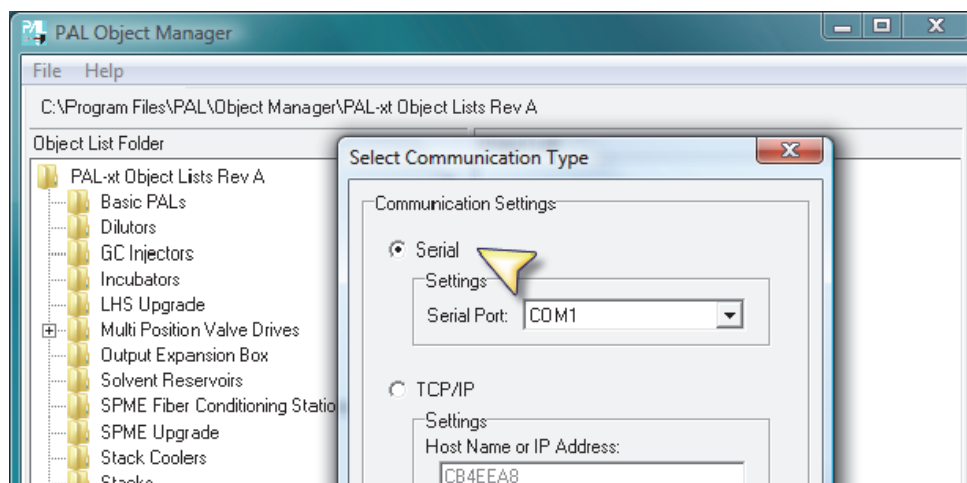
### ❖ To set up the communication mode

1. Choose **File > Options > Communication** or **Object List folder** to open the Select Communication Type dialog box.



2. Select the **Serial** option (Figure 99).

**Figure 99.** Select Serial communication



3. From the Serial Port list, select the COM $n$  port to use for communication.
4. Click **OK** to set the serial communication.

If no error message appears, the communication with the Accela Open Autosampler has been established. The Status Bar displays this message:

Connected on COM $n$

where  $n$  is the number of the COM port.

## Troubleshooting the Serial Communication Mode

PAL Object Manager software version 2.1.0. can address only COM port numbers 1 to 4. The ports must be consecutive. If you connect to one port and do not establish serial communications, connect to another port. Accela Open Autosampler

This limitation has been eliminated with PAL Object Manager software version 2.2.x.

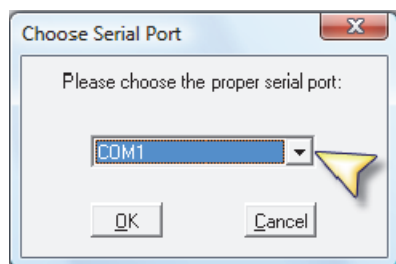


If you cannot establish serial communication with the PAL Object Manager software, check the following questions:

- Has the original RS232 cable been used, as provided with Accela Open Autosampler? (The PAL RS232 cable has a crossed wiring connection.)  
P/N 00950-01-00332 (3 m length) or P/N 00950-01-00333 (6 m length).
- Is the COM port on the computer available and correctly addressed?
- Are default settings verified?
- Is the COM port setting in the PAL Object Manager software correct?

If not, the message shown in [Figure 100](#) appears.

**Figure 100.** Wrong COM Port selected



These types of problems are seen mostly when a USB/Serial port is used to overcome the lack of a serial interface, as in today's laptop computers.

**Note** These two limitations surrounding the COM port setting are valid for the previous Object Manager software, Version 2.1.0. With the release of the new Version 2.2.x these problems are eliminated. The software scans for the available COM ports, the error message as shown in [Figure 100](#) is no longer displayed.

If no RS232 (serial) port is available on the computer, use a USB/Serial converter.

## Using PAL Object Manager to Change Your System Configuration

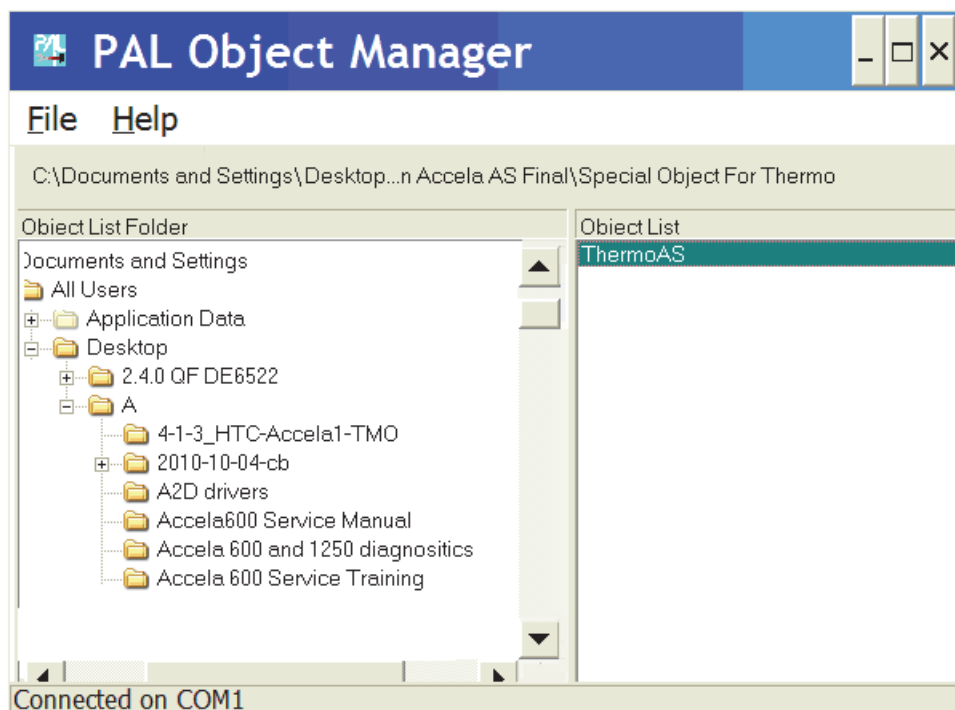
This section describes how to change your system configuration using the PAL Object Manager.

- [Starting the PAL Object Manager Application](#)
- [Using the Object Manager Application](#)

## Starting the PAL Object Manager Application

❖ **To start the PAL Object Manager application**

1. From the desktop, click the **PAL Object Manager** icon (  ).



The application opens and displays the various user buttons. The main part of the application shows the Object List folder pane with the current Object Lists folder.

All objects must correspond with the proper version of the firmware as identified by Thermo Scientific. For example, for objects using 4.x.x firmware, valid revisions are A, B, or C. For objects using a previous version of the software, use revision K.

The status bar displays the previously used communication mode.

2. To see the version information, choose **Help > About**.

The software displays the Object Manager software version.



## Using the Object Manager Application

The main purpose of the PAL Object Manager software is to load PAL firmware object lists on a system that has added a hardware module. To add an object list, follow this procedure, which describes adding a syringe object to an existing Accela Open Autosampler.

You can only add certain objects to your system. Be sure you know the impact of an object before adding it to the system.

Ensure that the Accela Open Autosampler is set up properly, the PAL Object Manager software is installed, and the communication is established as shown in “[Setting Up the Communication Mode](#)” on page 182.

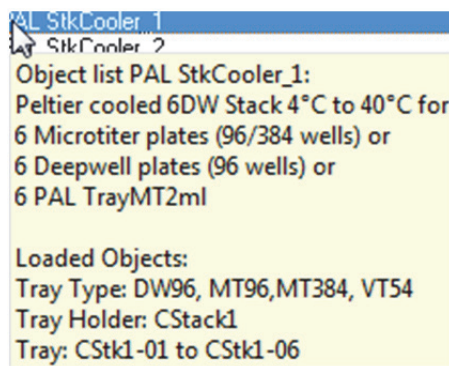
❖ **To add an object list to an existing Accela Open Autosampler**

1. In the Object List folder pane, click the **Syringes** folder.
2. Double-click the subfolder containing syringe firmware to open up the list of various available LC syringes.
3. In the Object List, select a specific syringe to highlight it.
4. Hold the cursor on the object list item to reveal information about the object.

This information provides the title and an explanation of the selected object.

The Loaded Objects area lists the objects that are loaded with one action. In this example it is a single object, but with other examples, such as a stack cooler, the situation is more complex ([Figure 101](#)). Four tray types, a tray holder (CStack1) and six trays (01-06) are sent to the Accela Open Autosampler in one package.

**Figure 101.** Object list stack cooler 1 example



Review the information to make sure that the correct object list has been selected and to understand what is included in this list. For example, you need to know whether all necessary tray types for the application are included or if you must specifically select another type and send it to the Accela Open Autosampler as a second task.

5. Click **Send to** Accela Open Autosampler to read the Accela Open Autosampler configuration and send the selected objects to the Accela Open Autosampler.

The configuration file is temporarily stored in the computer memory.

After the file has been sent successfully to the Accela Open Autosampler, the message “xx objects received” is displayed in the status bar, providing feedback from the Accela Open Autosampler to the Object Manager application.

In the example using the LC syringe, 25  $\mu\text{L}$ , the Object Manager software might send two objects in a single file. In this case, the first object contains the geometric definition of the syringe itself and the second object specifies which plunger motor driver is assigned to this particular syringe.

The same concept is used to send the desired objects to the Accela Open Autosampler for all other object lists.

The Object Manager always lists the object being sent in the last line of the PAL firmware class, visible from the local terminal. The Object Manager software can quickly check whether the PAL firmware has received the new object.

In some cases—seen in particular in the PAL but not in the PAL-xt System—you might receive an error message: Memory full. To resolve this problem, you must delete as many objects as possible from the Accela Open Autosampler that you are not using. You can link a syringe to a method and a job. First delete the job, then the method, and finally the syringe object. See “[To delete objects](#)” on [page 186](#) for information about deleting objects.

Confirm your understanding of the Object List parameters and descriptions to make sure that the correct, desired objects are sent to the PAL firmware. In certain cases where complexity is greater—for example, multi-position valve drives—the guideline provides the correct order of steps in the process to provide reliable functioning of the Accela Open Autosampler.

#### ❖ **To delete objects**

Use the Accela Open Autosampler handheld device to delete objects.

In the handheld device, objects are displayed as a tree. Identify the object and delete it according to instructions in “[Operating the Accela Open Autosampler](#)” on [page 55](#).

To delete multi-object items, delete the lowest items first and then remove higher objects until you delete the top-level object.

## Column Switching Option

This appendix describes how to install and use the column selector option for the Accela Open Autosampler.

### Contents

- [Installing the Column Switching Valves](#)
- [Installing the Software Tools](#)
- [Backing Up the System](#)
- [Installing the Column Switch Option](#)
- [Verifying the Valve Configuration](#)
- [Installing the Macros and Methods](#)
- [Creating a Method in Xcalibur](#)

## Installing the Column Switching Valves

To install the 6-port and 10-port valves for the column selector option, follow these procedures:

- [Installing the 6-Port Valve](#)
- [Installing the 10-Port Valve](#)

## Installing the 6-Port Valve

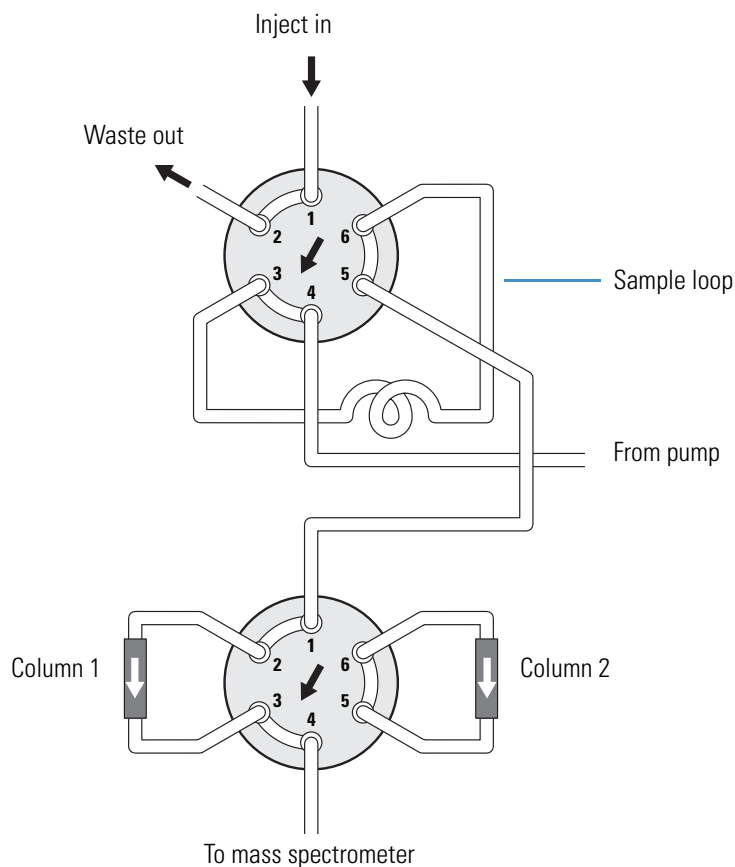
### ❖ To install the 6-port valve

1. Install the valve drive onto the valve bracket that is attached to the *x*-axis rail.

**Note** To minimize gradient delay volume, install the 6-port valve as close as possible to the injection valve as possible. For best results, you use the middle position of the 3-valve holder stack.

2. Connect the MODBUS A connector on the valve drive to the MODBUS connector on the back of the *x*-axis main rail using the supplied cable assembly. If several serial valve drives are installed, connect each valve drive in serial.
3. Install the 6-port valve into the valve drive and secure in place using a T-20 Torx driver, with port 1 facing up.
4. Install all tubing and columns using the tubing, fittings, and ferrules supplied in the kit (Figure 102).

**Figure 102.** Two column switching using 6-port valve



## Installing the 10-Port Valve

This section describes how to install the 10-port valve for the Column Switch Option.

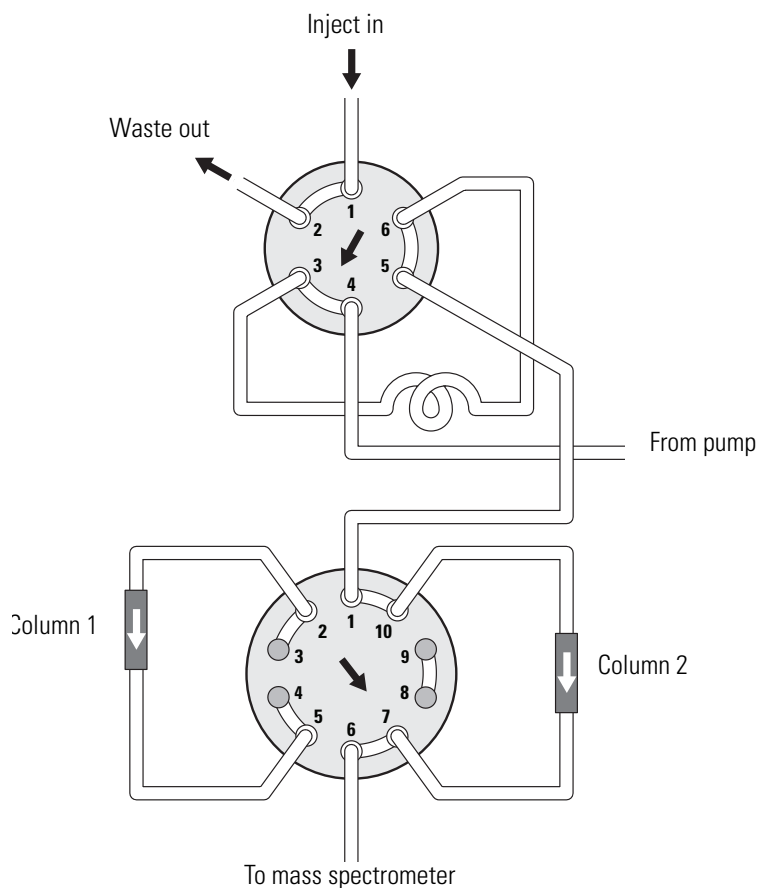
### ❖ To install the 10-port valve

1. Install the valve drive onto the valve bracket that is attached to the *x*-axis rail.

**Note** The location of the installed 10-port valve should be as close to the injection valve as possible to minimize gradient delay volume. It is recommended that you use the middle position of the 3-valve holder stack.

2. Connect the MODBUS A connector on the valve drive to the MODBUS connector on the back of the *x*-axis main rail using the supplied cable assembly.
3. Install the 10-port valve into the valve drive and secure in place using a T-20 Torx driver, with port 1 facing up.
4. Install all tubing and columns using the tubing, fittings, and ferrules supplied in the kit (Figure 103).

**Figure 103.** Two column switching using 10-port valve

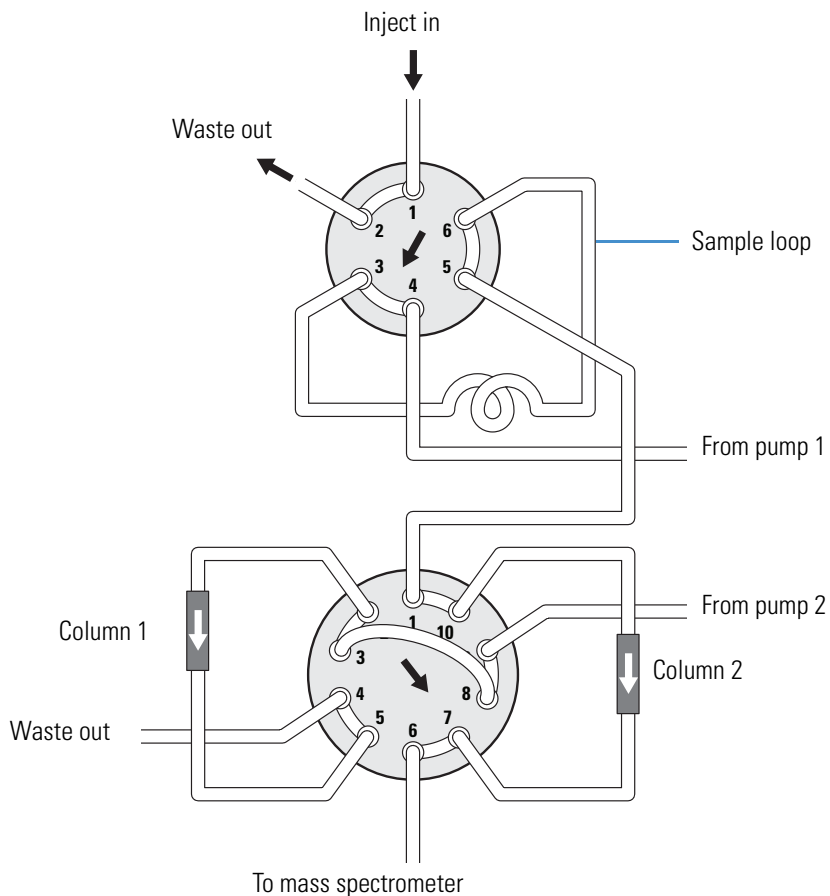


Note: Plug ports 3, 4, 8, and 9 in the 10-port valve.

**IMPORTANT** For switching between the two columns, make sure that ports 3, 4, 8 and 9 are plugged to prevent contamination from entering these ports.

For plumbing the system as shown in [Figure 104](#) (for column conditioning), you must install two Accela pumps. Refer to the *EQuan Getting Connected Guide* for the installation procedures.

**Figure 104.** Two column conditioning using 10-port valve (Note: Requires two pumps)





## Installing the Software Tools

This section describes the procedure to install the software tools for the column switch option.

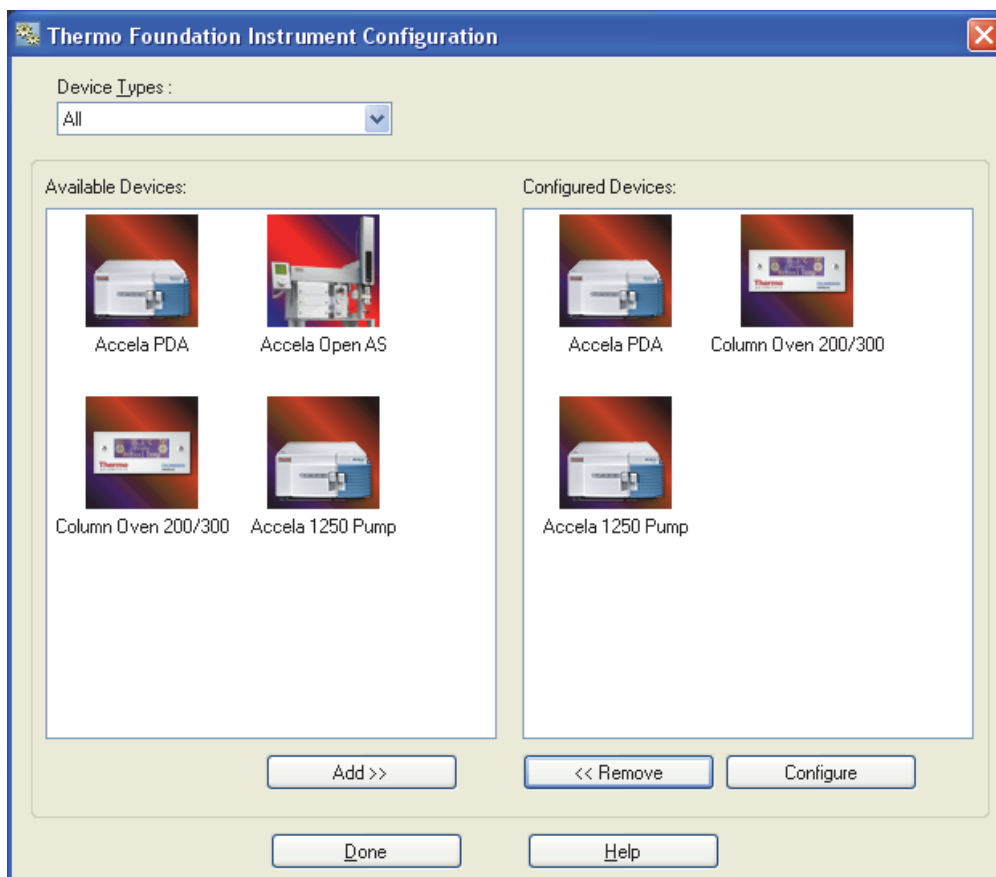
❖ **To install the software tools**

1. Install the PAL Loader as described in “[Installation of PAL Loader Software](#)” on [page 156](#). Make sure to use PAL Loader version 2.1.0 or later.
2. Install the Object Manager as described in “[Installing PAL Object Manager](#)” on [page 180](#). Make sure to use Object Manager 2.2.1.

**Note** The PAL Loader and Object Manager are included on the Accela Open Autosampler Controls software CD that comes with your Accela Open Autosampler.

**Note** If Xcalibur 2.1 or later is installed, remove “Accela Open AS” from the Configured Devices list in the Instrument Configuration window ([Figure 105](#)).

**Figure 105.** Thermo Foundation Instrument Configuration window



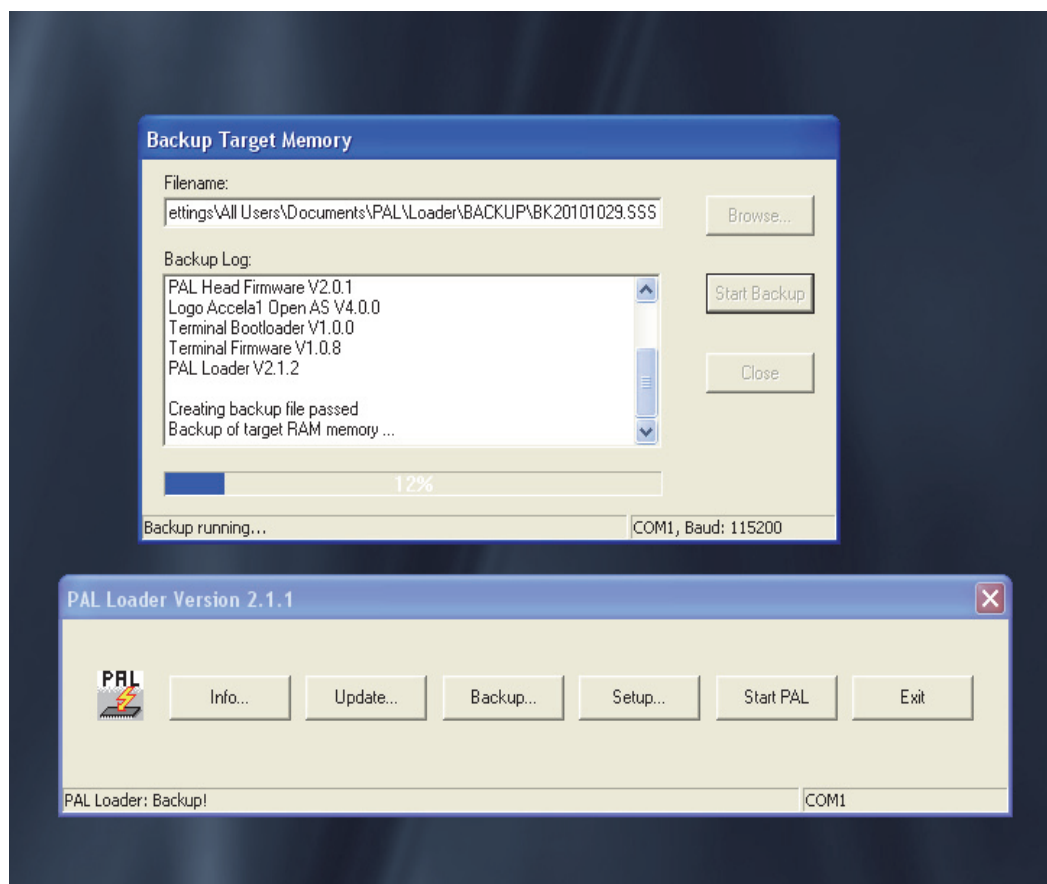
## Backing Up the System

This section describes how to create a backup file for the system.

Before you install the column switch option or make any changes to the autosampler, always create a backup file using PAL Loader 2.1.0 or later so that you can restore the system in case of a crash or error.

Create the backup file as shown in [Figure 106](#).

**Figure 106.** Backup file creation



**Tip** Thermo Fisher Scientific recommends that you create a new backup file after the system has worked with the column switching objects.

## Installing the Column Switch Option

Use one of the following two methods to install the column switch option:

- [Installing a New Backup File to the Autosampler](#)
- [Installing Special Objects to the Autosampler](#)

**Note** If you install the new backup file, you must teach all positions again. If you choose to install the special objects instead of the new backup file, you don't need to teach all positions again, but you must manually install all new objects. The new backup file installs these new objects for you.

### Installing a New Backup File to the Autosampler

**IMPORTANT** If you use this method, you must teach all positions again.

To install a new backup file, follow these procedures:

1. [Reading Out and Recording the Serial Number](#)
2. [Installing Option from the Backup File](#)
3. [Re-entering the Serial Number](#)

#### Reading Out and Recording the Serial Number

Before installing the new backup file, you must read out and record the existing serial number on the control terminal (hand-held controller) of the Accela Open Autosampler.

❖ **To read out the serial number**

1. Choose **Menu > F3-Setup > Ident > System SNo.**
2. Read out the serial number and record it for later use.

#### Installing Option from the Backup File

❖ **To install option from the backup file**

1. On the CD that comes with the column switch kit, locate the following files:

For Firmware 4.1.2, go to *drive:\Backup\Firmware 4.1.2\*

Valve option	Open AS Config 1 (3 drawer)	Open AS Config 2 (6 drawer)
6-port valve option	4-1-2_HTC-Accela1-TM O-ad-6PortS_B.SSS	4-1-2_HTC-Accela2-TM O-ad-6PortS_B.SSS
10-port valve option	4-1-2_HTC-Accela1-TM O-ad-10PortS_B.SSS	4-1-2_HTC-Accela2-TM O-ad-10PortS_B.SSS

For Firmware 4.1.3, go to *drive:\Backup\Firmware 4.1.3\*

Valve option	Open AS Config 1 (3 drawer)	Open AS Config 2 (6 drawer)
6-port valve option	4-1-3_HTC-Accela1-TM O-ad-6PortS_C.SSS	4-1-3_HTC-Accela2-TM O-ad-6PortS_C.SSS
10-port valve option	4-1-3_HTC-Accela1-TM O-ad-10PortS_C.SSS	4-1-3_HTC-Accela2-TM O-ad-10PortS_C.SSS

2. Copy and save the files in the tables above to the folder: C:\Thermo\Instruments\LC Devices\AccelaOpenAS on your computer.
3. Check for the correct configuration. Config 1 is for use with the 3-drawer cooled stack. Config 2 is for use with the 6-drawer cooled stack.
4. Update the autosampler with PAL loader 2.1.1 or later with the correct SSS file.
5. Teach all position after uploading the new SSS file as described in the *Accela Open Autosampler Hardware Manual*.

## Re-entering the Serial Number

After installing option from the backup files, re-enter the serial number using the control terminal (hand-held controller).

### ❖ To re-enter the serial number

1. Choose **Menu > F3-Setup > Ident > System SNo.**
2. Re-enter the system serial number.

## Installing Special Objects to the Autosampler

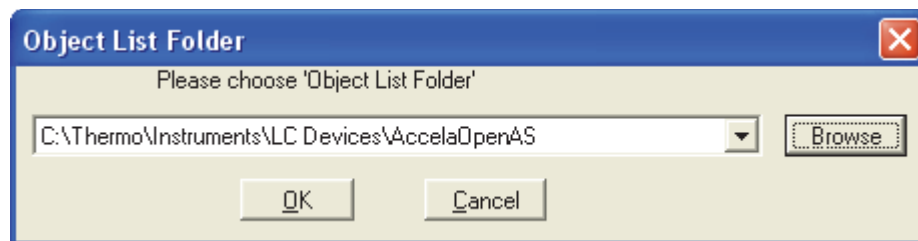
**IMPORTANT** Follow these instructions only if you choose to install the special objects and do not install the new backup file to the autosampler using the PAL Loader.

### ❖ To install special objects to the autosampler

1. On the CD that comes with the column switch kit, locate the following files:
  - For Firmware 4.1.2, go to:  
*drive:\Objects\Firmware 4.1.2\PAL Acc-C72VX6port\_B.pol*  
*drive:\Objects\Firmware 4.1.2\PAL Acc-C72VX10port\_B.pol*
  - For Firmware 4.1.3, go to:  
*drive:\Objects\Firmware 4.1.3\PAL Acc-C72VX6port\_C.pol*  
*drive:\Objects\Firmware 4.1.3\PAL Acc-C72VX10port\_C.pol*

2. Copy and save the files above in the following folder on you computer:  
C:\Thermo\Instruments\LC Devices\AccelaOpenAS
3. Check for the correct firmware version. B indicates object files for firmware 4.1.2.  
C indicates object files for firmware 4.1.3 and later.
4. Start the Object Manager.
5. Browse and set the object list folder as shown in [Figure 107](#).

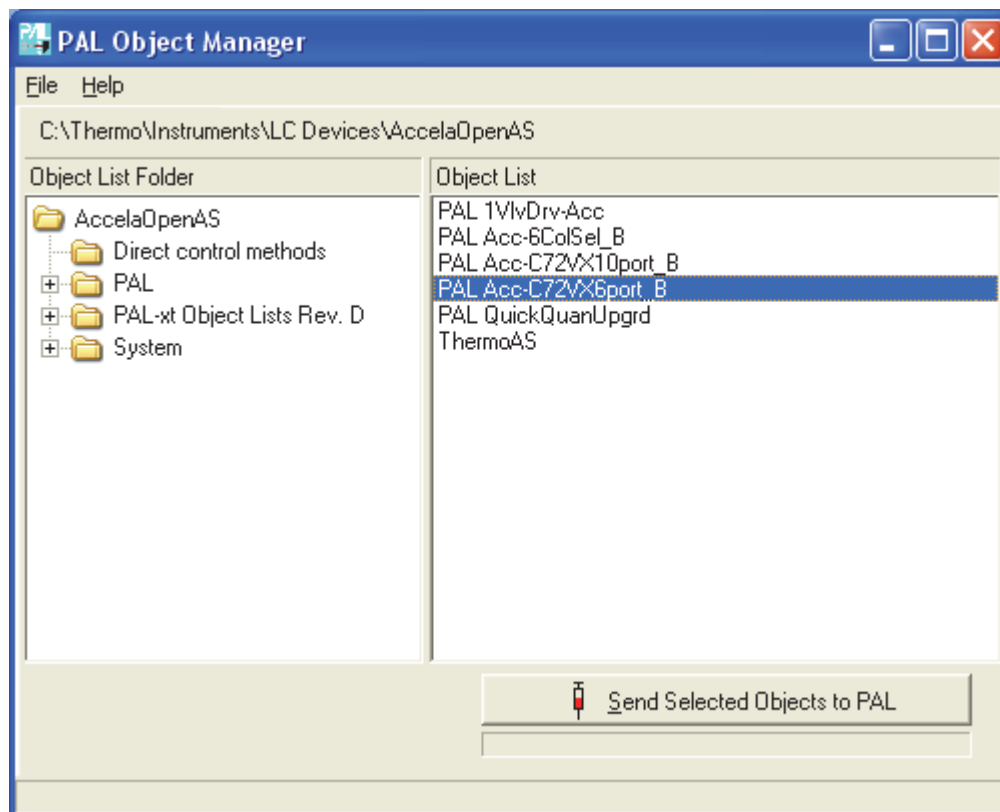
**Figure 107.** Object list folder setting



6. Select **PAL Acc-C72VX6port\_X** or **PAL Acc-C72VX10port\_X**, and click **Send Selected Objects to PAL** ([Figure 108](#)).

**Note** X=B for firmware 4.1.2; X=C for firmware 4.1.3.

**Figure 108.** PAL Object Manager dialog box



7. Restart the Accela Open Autosampler by turning the power switch off and then on.
8. If the “Ser Valve MV2 Not Detected” error message displays, do one of the following:
  - Check if the MODBUS connectors on the back of the Accela Open Autosampler and the valve drive are secure; then check if the additional valve drive is detected using the control terminal (hand-held controller) as described below.
    - i. On the top level of the control terminal screen, press F1 for Menu, and then choose **Utilities > Injector > LC Vlv4**.
    - ii. Press F2 (Act Valve) to rotate the valve, noting that the red LED illuminates.
    - iii. Press F2 (Deact Valve) to rotate the valve back to its original position (the LED extinguishes).
    - iv. Press F4 to go back to the Home screen.

–or–

- Check if the ID number of the connected valve is correct. If not, set the correct ID number (located on the back of the valve drive) using the control terminal (hand-held controller) as described below. Be sure that only one valve drive is connected to the MODBUS connector.
  - i. On the top level of the control terminal screen, press F1 for Menu, and then choose **Setup > Objects > Serial valves > Set VlvID**.
  - ii. Press F4 to go back to the Home screen.

## Verifying the Valve Configuration

**Note** Regardless of which of the two methods you use to install the column switch, you must verify the valve configuration as described in this section.

After installing the column switch option, use the control terminal (hand-held controller) to verify the valve configuration settings as shown in [Table 50](#). Make sure the settings shown on the control terminal screen match the parameters in [Table 50](#).

### ❖ To verify the valve configuration using the control terminal

Choose **Menu > F3-Setup > Objects > Serial Valves**.

**Note** The LED on the front of the valve drive illuminates red when it is in the On position.

**Table 50.** Valve configuration parameters

Setting	6 port valve	10 port valve
Max Position	6	10
Off Position	1	1
On Position	2	2
Mode	HSpeed	HTorque

## Installing the Macros and Methods

### ❖ To install the macros and methods

1. On the CD that comes with the column switch kit, locate the following files and copy them to the appropriate folders:

Find these files:

*drive:\Macros and Cycles\Xcalibur\Macros\Standard Inj Accela Open Col 1.pma*

*drive:\Macros and Cycles\Xcalibur\Macros\Standard Inj Accela Open Col 2.pma*

Copy these two files to this folder on your computer:

C:\Thermo\Instruments\LC Devices\AccelaOpenAS\PAL\Macros

Find these files:

*drive:\Macros and Cycles\Xcalibur\Methods\Standard Injection Accela Open Col1.pme*

*drive:\Macros and Cycles\Xcalibur\Methods\ Standard Injection Accela Open Col2.pme*

Copy these two files to this folder on your computer:

C:\Thermo\Instruments\LC Devices\AccelaOpenAS\PAL\Methods

Find these files:

*drive:\Macros and Cycles\Xcalibur\Direct Control Methods\Direct Control - Switch Column Col 1.meth*

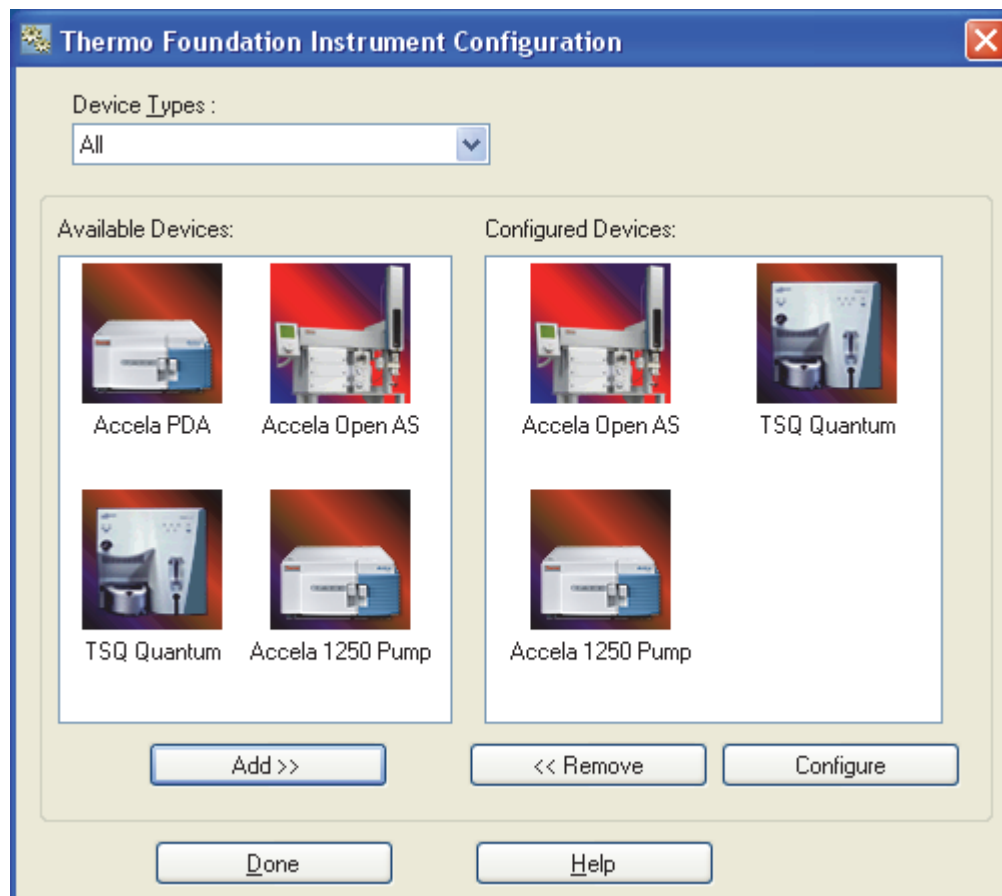
*drive:\Macros and Cycles\Xcalibur\Direct Control Methods\Direct Control - Switch Column Col 2.meth*

Copy these two files to this folder on your computer:

C:\Thermo\Instruments\LC Devices\AccelaOpenAS\PAL\DirectControlMethods

2. Add the Accela Open Autosampler back into the Configured Devices in the Instrument Configuration as shown in [Figure 109](#).

**Figure 109.** Thermo Foundation Instrument Configuration



## Creating a Method in Xcalibur

After installing the valves, the software tools, the column switch option, and the macros and methods, create a method in the Xcalibur data system to test the column selector option.

### ❖ To create a method the Xcalibur data system

1. From the Xcalibur Home Page, click the **Instrument Setup** icon, .

The Instrument Setup window opens with icons for the configured instruments in the viewbar on the left side of the window.

2. In the viewbar, click the **Accela Open AS** icon, .

The Accela Open AS view opens.

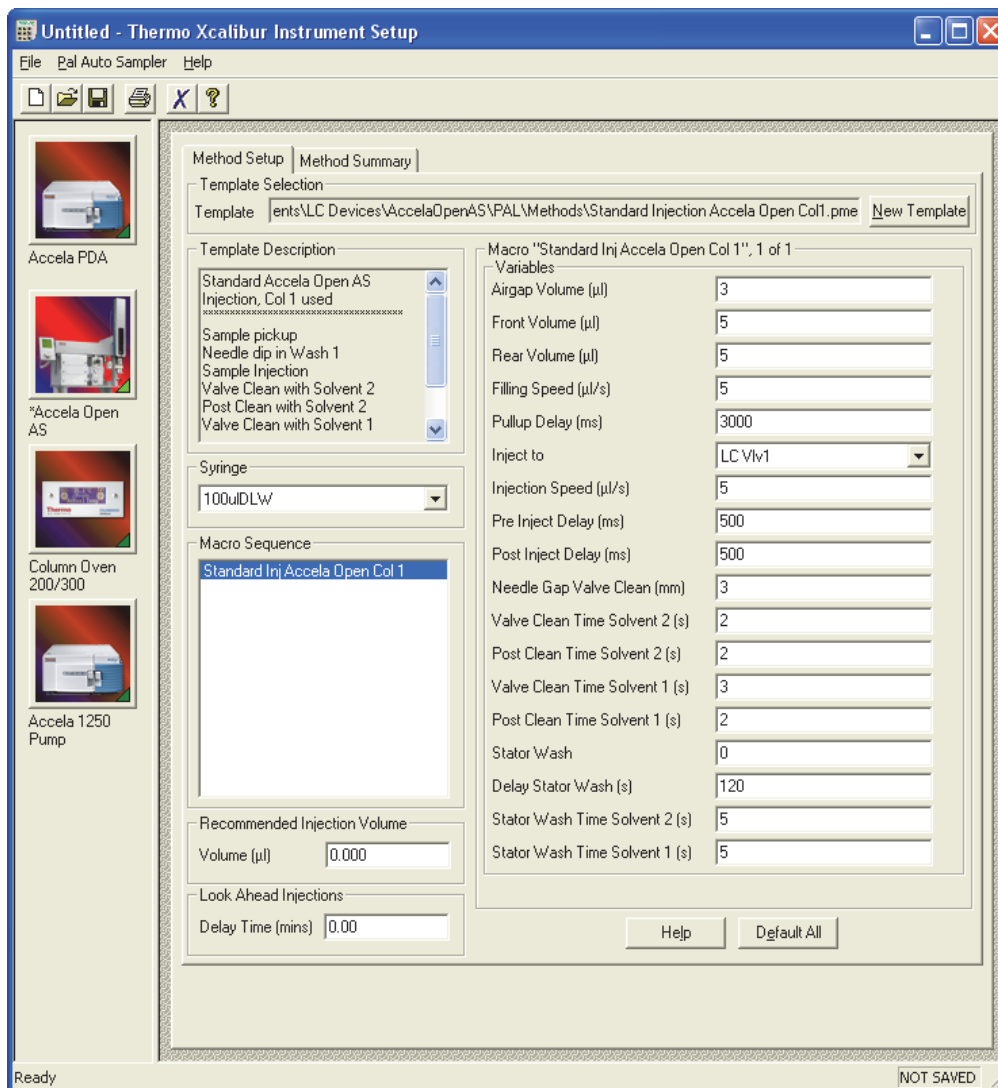
3. Click the **Method Setup** tab.



4. Create the method as shown in [Figure 110](#). Make sure you do the following:

- Use the “Standard Injection Accela Open Col1” method for injection to column 1.
- Use the “Standard Injection Accela Open Col2” method for injection to column 2.

**Figure 110.** Thermo Xcalibur Instrument Setup window



**IMPORTANT** As a precaution, do the following after creating the method:

- Check the flow paths for column 1 and column 2 after installation by testing the flow paths with liquid.
- Create labels and place them on the fluidic connections to identify entry ports for column 1 and column 2.
- Back up the system again as shown in “[Backing Up the System](#)” on [page 192](#).



## Changing the DLW Holding Loop

With the DLW-2 option, a new DLW holding loop design has been introduced. The DLW-2 holding loop is a high performance loop, which is compatible with the previously used FEP loop with a separate needle and flow diverter. The DLW-2 holding loop gives better performance in regard to carryover and low volume injection.

Table 51 shows the differences between the DLW and DLW-2 holding loops.

**Table 51.** Comparison of the two holding loop designs (Sheet 1 of 2)

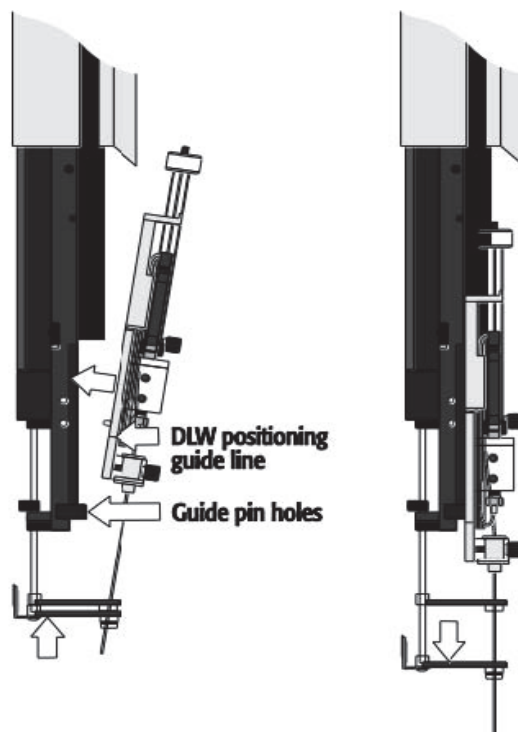
Part or Parameter	DLW holding loop	DLW-2 holding loop
<b>Holding loop design</b>	P/N 00950-01-00317	<b>P/N 00950-01-00360</b>
Material:	FEP tubing	Loop stainless steel, passivated
OD:	1/16 in. (1.58 mm)	0.72 mm (Gauge 22)
ID:	0.50 mm	0.41 mm
Length:	550 mm	855 mm
Volume:	108 µL	118 µL
<b>Syringe needle</b>	P/N 00950-01-00328	Not applicable, integrated in DLW-2 holding loop
Length:	51 mm	
Gauge:	22	
Material:	Stainless steel	
<b>DLW flow diverter</b>	P/N 00950-01-00319	Not applicable, integrated in DLW-2 holding loop
Length:	20.5 mm	
OD:	0.72 mm (Gauge 22)	
ID:	0.41 mm	
Material:	Stainless steel	

**Table 51.** Comparison of the two holding loop designs (Sheet 2 of 2)

Part or Parameter	DLW holding loop	DLW-2 holding loop
<b>DLW internal volume</b>		
Total delay volume:	205 $\mu\text{L}$ + injection loop volume	208 $\mu\text{L}$ + injection loop
	Manifold: 90 $\mu\text{L}$	Manifold: 90 $\mu\text{L}$
	Holding loop: 108 $\mu\text{L}$	Holding loop: 118 $\mu\text{L}$
	Syringe needle: 6.7 $\mu\text{L}$	
	Installed injection loop	Installed injection loop

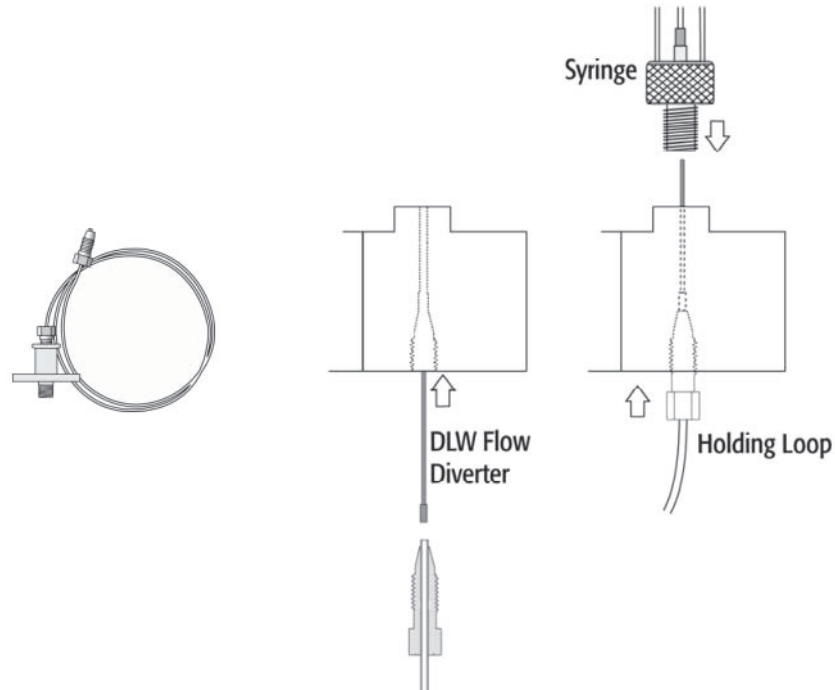
❖ **To change the DLW-2 holding loop**

1. Turn off the PAL-xt instrument.
2. Pull down the injection unit slider as far as possible in order to gain access to DLW syringe and holder assembly.
3. Remove the DLW syringe from the DLW-2 Syringe Holder Assembly.
4. Unscrew the knurled screw on the holder and remove the DLW-2 Syringe Holder Assembly from the Injection Unit. Thereby hold up the lower needle guide to prevent the needle from bending (Figure 111).

**Figure 111.** Remove the DLW-2 syringe holder assembly

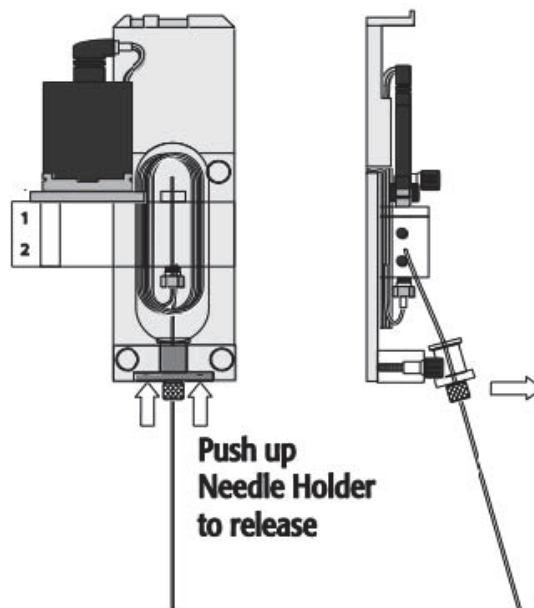
- If there is a FEP holding loop installed, loosen the retaining nut holding the needle to the needle adapter (Figure 112).

**Figure 112.** Loosen the retaining nut for the FEP holding loop



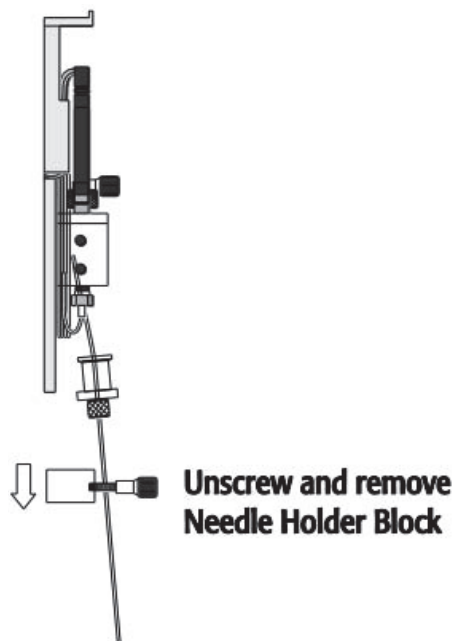
If there is a DLW-2 holding loop installed, push up the DLW needle holder assembly and click it out of the DLW Needle Adapter Block (Figure 113).

**Figure 113.** Push up the needle holder to release



6. Unscrew and remove the Needle Holder Block (Figure 114).

**Figure 114.** Unscrew and remove the Needle Holder Block

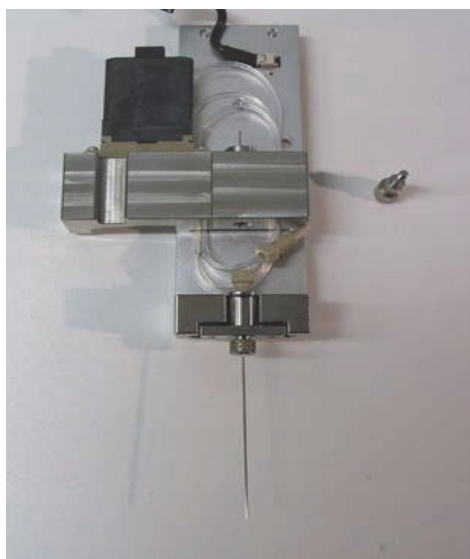
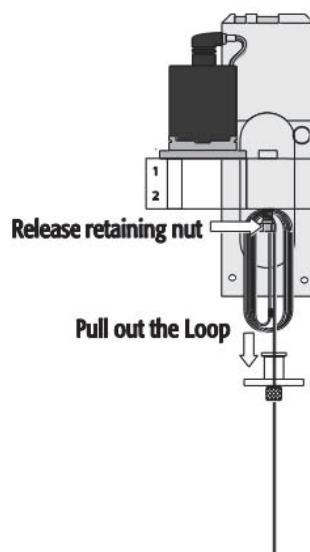


7. Unscrew the retaining nut that tightens the holding loop to DLW Manifold (both designs). Pull out the DLW holding loop (Figure 115).

**Figure 115.** Release the retaining nut and pull out the loop

DLW-2 needle loop

FEP (DLW legacy) needle loop



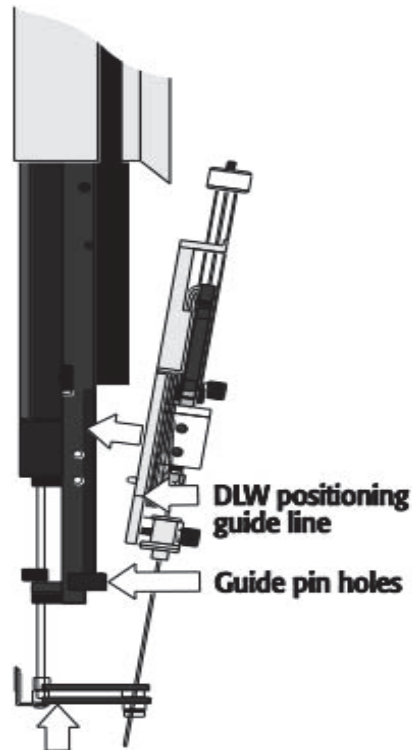
If there is an FEP holding loop installed, remember to remove now the flow diverter from the DLW Manifold (Figure 116). Removing the flow diverter is not necessary if the DLW-2 holding loop is installed because the flow diverter is part of the holding loop.

**Figure 116.** Remove the flow diverter



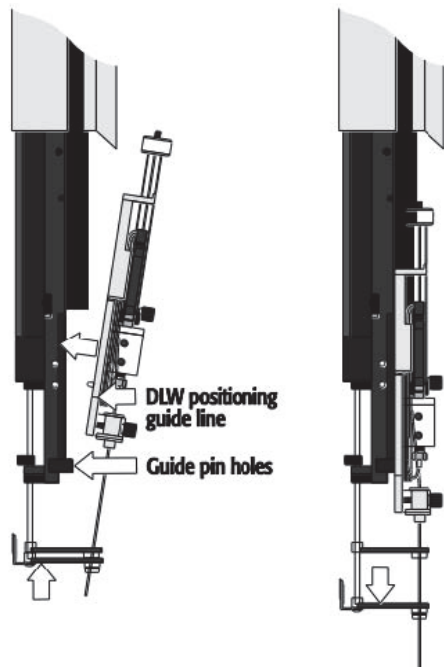
8. Insert the DLW-2 holding loop to the DLW Manifold and tight the ferrule.
9. Fix the DLW-2 holding loop with the slotted needle adapter to the DLW-2 Syringe Holder Assembly. Make sure the opening of the needle adapter is facing back.
10. Fix the DLW-2 holding loop to the needle adapter with the retaining nut.
11. Insert the DLW-2 syringe adapter without installed syringe (Figure 117).

**Figure 117.** Install the DLW-2 syringe adapter



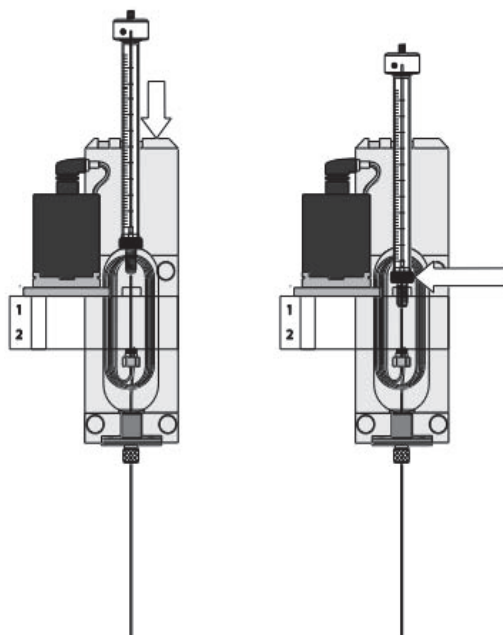
12. After the DLW-2 syringe holder is mounted on the Z-Unit, tighten the knurled screw to fix the holder assembly on the syringe slider (Figure 118).

**Figure 118.** Fix the holder assembly on the syringe slider



13. Insert the DLW syringe and tighten it (Figure 119).

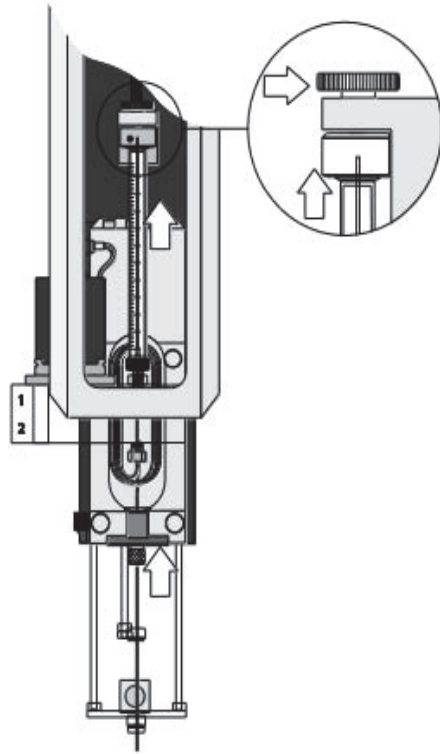
**Figure 119.** Insert the DLW syringe





14. Connect the DLW syringe plunger holder (Figure 120).

**Figure 120.** Connect the DLW syringe plunger holder







# Index

## A

Accela Open Autosampler  
assembling 14  
components 1  
electrical connections 36  
error messages 85  
external connectors 145  
firmware overview 136  
installing 11  
maintenance 81  
naming convention 116  
object positions 40  
operating 55  
operating and environmental requirements 9  
physical specifications 8  
replacement parts 101  
sound pressure level 9  
specifications 6  
system limitations 152  
unpacking components 12  
using control terminal 55

## B

backing up PAL firmware, PAL Loader 166

## C

column switch option 187  
backing up system 192  
installing 193  
installing macros and methods 197  
installing software tools 191  
installing the 10-port valve 189  
installing the 6-port valve 188  
verifying valve configuration 196  
communication mode, setting up for PAL Object Manager software 182

compliance  
FCC iii  
regulatory iii  
compliance, WEEE v  
contacting us xvii  
control terminal  
function keys 57  
illustration 56  
menu screens 56  
using 55  
Cycle 114

## D

DLW  
Actuator/Solenoid 27  
characterizing features 3  
Flow Diverter 12  
Holding Loop 12–13  
manifold 27  
needle guide length tool 12, 14  
overview 3  
Plunger Holder 12–13  
Pump Module 12–13  
replacement 101  
spare parts ordering information 103  
Syringe 12, 29  
Syringe Holder assembly 12–13  
Syringe Needle 12–13  
Tubing Kit 12–13  
DLW pump module, replacing 102  
DLW replacing parts  
syringe 102  
syringe holder assembly 102  
wash station 102  
DLW Syringe Holder Assembly, installing 27  
DLW Wash Station  
Injection Valve plumbing and connection 35  
installing 25

## Index: E

DLW-2  
  components 13  
  definition 4  
  spare parts ordering information 108  
documentation  
  accessing xvi  
  additional xv  
  online xvii  
  survey xvii

Dynamic Load and Wash (DLW) 3

## E

electromagnetic compatibility iii  
EMC compliance iii  
Extended User Mode, accessing 117  
external connectors, Accela Open Autosampler 145

## F

Fast Wash Station for two different solvents 2  
FCC compliance iii  
firmware  
  error codes 140  
  overview 136  
flow path, injection valve 50

## I

info functions 68  
  free objects/free items 70  
  hardware 69  
  maintenance 69  
  software 69  
Injection Unit 2  
  installing 18  
  ribbon cable 19  
injection unit 12  
injection valve  
  flow path 50  
  Valco W-Type 6-Port 51  
introduction, Object Manager Application 184

## J

Job 114  
Job queue 114  
job queue 59  
  building and starting 60  
  canceling 61

## K

keypad terminal 2, 12  
keypad terminal, installing 21

## L

LC injection valve, installing 22  
logfile 68

## M

Method 114  
Method parameters 114  
methods 57  
  creating 57  
  deleting 59  
  editing and viewing 58  
Module 114

## O

Object class 114  
Object Item 114  
Object list 115  
Object Manager 115  
Object Manager Software, installing 180  
object positions, descriptions 41  
Objects 114  
Open Autosampler components 1  
ordering information  
  DLW spare parts 103  
  DLW-2 spare parts 108

## P

PAL Firmware, updating 168  
PAL Loader 155  
  backing up PAL Firmware 166  
  dedicated shortcut for serial communication 159  
  installing 156  
  multiple shortcuts assignment 158  
  new features 156  
  operating 161  
  reset PAL-xt to serial communication 165  
  serial communication 164  
  setting up COM ports 162  
  starting 161  
PAL Object Lists folder  
  installing for PAL Object Manager software 181  
  PAL Object Manager Software 184  
PAL Object Manager software  
  backing up your system configuration 179  
  installing 180

- port settings
  - choose COM port for PAL Object Manager software 182
  - troubleshooting for PAL Object Manager software 183
- power supply, installing 22
- pump holder, mounting 23
- pumps, self-priming 3

## R

- reference points
  - injection valve 48
  - Wash Station 46
  - waste 47
- regulatory compliance iii
- resetting PAL-xt to serial communication, PAL Loader 165

## S

- safety guard 2, 12
- safety standards iii
- self-priming micro pumps 3
- serial communication, PAL Loader 164
- setting up COM parts, PAL Loader 162
- setting up LAN communication, PAL Loader 165
- setup functions
  - objects 72
  - selecting 71
  - sounds 71
  - time 71
- Stack 115
- Stack Cooler
  - condensation build-up 78
  - temperature alarm 77
  - temperature control 76
  - temperature setting 75
  - temperature stability 77
- stand-alone supports 12
- standard chromatographic tests 96
- standard liquid syringe holders 27
- standard stand-alone supports 2
- starting, PAL Object Manager Application 184
- survey link xvii
- syringe kit 2, 12
- system components, Accela Open Autosampler 2
- system requirements, PAL Object Manager software 180

## T

- Torx driver 19
- Tray 115

- tray
  - holder 43
  - position 43
  - rows and columns 43
- Tray holder 115
- Tray type
  - parameter 115
  - pattern explanation 131
- troubleshooting
  - Accela Open Autosampler 85
  - PAL Loader 171
  - special considerations 90

## U

- updating PAL firmware, PAL Loader 168
- utility functions
  - description 61
  - dilutors 66
  - injector 64
  - syringe 62
  - tools 67
  - tray 63
  - vial 66
  - wash station 65

## V

- Valve Needle Guide Assembly, installing 34
- valve penetration value, defining 48

## W

- Wash Station assembly 12
- Wash Station reference point 46
- WEEE compliance v

## X

- $x$ -,  $y$ -axes assembly 2

## Z

- $z$  axis 2

