

EASY-nLC II

Version 2.8

User Guide

60053-97000 Revision A September 2010

DOCUMENTATION
SURVEY

© 2010 Thermo Fisher Scientific Inc. All rights reserved.

EASY-nLC, EASY-Column, AFC, and IFC are trademarks, and Microtiter is a registered trademark of Thermo Fisher Scientific Inc.

PEEKsil is a trademark of SGE International Pty Ltd Corp.

This product uses FreeRTOS from www.freertos.org (available upon request). FreeRTOS is a trademark of Real Time Engineers Ltd.

The following are registered trademarks in the United States: Advion and RePlay are registered trademarks of Advion BioSystems, Inc. Duran is a registered trademark of Schott AG. PhotoMOS is a registered trademark of Panasonic Electric Works, Co., Ltd. VICI and Valco are registered trademarks of Valco Instruments Co., Inc.

Microsoft and Windows are registered trademarks of the Microsoft Corporation in the United States and other countries.

All other trademarks are the property of Thermo Fisher Scientific 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.

Thermo Fisher Scientific Inc. makes no representations that this document is complete, accurate or error-free and assumes no responsibility and will not be liable for any errors, omissions, damage or loss that might result from any use of this document, even if the information in the document is followed properly.

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, September 2010

Software version: EASY-nLC II version 2.8

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

Contents

	Preface	vii
	Related Documentation	viii
	Safety and Special Notices	viii
	Safety Practices	viii
	Special Notices	viii
	Contacting Us	ix
	Declaration of Conformity	xiv
	U.S. Safety and EMC (Electromagnetic Compliance) Standards	xiv
	Canadian Safety and EMC (Electromagnetic Compliance) Standards	xv
	European Safety and EMC (Electromagnetic Compliance) Standards	xv
Chapter 1	Introduction	1
	Product Description	2
	System Components	2
	System Configuration for Two-Column Setup	3
	System Configuration for a One-Column Setup	4
	Instrument Operation	4
	Pump Flow Control	5
	Graphical Hardware Overview	7
Chapter 2	Installation	11
	Placement	11
	Power	11
	Temperature and Humidity	12
	Lifting Instruction	12
	External Connections	12
	Connecting to the Mass Spectrometer through Contact Closure	13
	Connecting to Remote Support	13
	Connecting to a Local Area Network for Data Exchange	14
	Attaching Mouse and Keyboard to the USB Connections	14
	Attaching Add-on Products through the P-Bus and RS-232	14
	Flow Parts and Lines	15
Chapter 3	Instrument Control Software	17
	Using the Touch-Sensitive Screen	18
	Starting the EASY-nLC II System	19

	User Interface Layout	20
	General Interaction Principles	21
	Application Menu Structure	21
	Home	22
	Batch Setup	23
	Method Editor	25
	Maintenance	27
	Configuration	29
	Logging In to the EASY-nLC II System.	32
	Closing Down the EASY-nLC II System.	33
Chapter 4	Configuring the EASY-nLC II System	35
	Setting Up the Mass Spectrometer Connection	35
	Network Connection.	36
	User Permissions	38
	Changing the Administrator Password.	39
	Creating New User Accounts.	40
	Configuring Column Setup, Loop Volume, Idle Flow, and IFC	40
Chapter 5	Preparation for Use	43
	Preparing Solvent Bottles and Waste Containers	43
	Solvent A and B Bottles.	44
	Autosampler Bottles	44
	Waste Container	45
	Executing Maintenance Scripts	46
	Purging and Flushing the Pumps.	47
	Installing Columns	50
	Preparing the Pre-Column.	51
	Equilibrating the Analytical Column.	52
Chapter 6	Running the First Sample	53
	Preparing a Well Plate	54
	Creating a Method	54
	Page 1/5	54
	Page 2/5	55
	Page 3/5	56
	Page 4/5	58
	Page 5/5	59
	Creating a Batch	60
	Starting Sample Acquisition	63
	Monitoring the Run	63
	Using the Overview Page in the Home Menu	64
	Using the Graphs Page in the Home Menu	65

	Stopping Sample Acquisition	66
	Editing the Running Batch	67
	Troubleshooting a Sample Run	70
Chapter 7	Autosampler Calibration	71
	Replacing the Autosampler Adapter Plate	72
	Managing Plate Formats	75
	Preparation and Basic Principles	77
	Calibrating Plates	78
	Calibrating the Wash/Waste Compartment.	81
Chapter 8	Troubleshooting.	87
	Running a Batch	88
	Autosampler	90
	Contact Closure	92
	Network Access	92
	Starting Up the System	93
	System Component States (Valves, Pumps, Autosampler, and Plate Cooler)	93
	Miscellaneous	94
Chapter 9	Remote Support	97
	Connecting the EASY-nLC II System to the Support Server	97
	Saving System Files on a USB Removable Storage Device	99
Chapter 10	Maintenance	103
	Daily Maintenance	103
	Quarterly Maintenance	104
	Yearly Maintenance	104
	Maintenance Scripts	105
	Prepare – Purge Solvent.	106
	Prepare – Flush Air	107
	Prepare – Precolumn Equilibration	108
	Prepare – Analytical Col Equilibration	109
	Prepare – Isocratic Flow	110
	Test – MS Connection	111
	Test – Sample Pickup	112
	Test – Leaks	113
	Test – Valve Check	114
	Test – Back Pressure	116
	Test – Autosampler Torque.	117
	Calibrate – Valve Tune	117
	Calibrate – Flow Sensors	117
	Calibrate – Reset Pressure Sensor.	118

Appendix A Specifications	119
Performance Specifications	119
Technical Specifications	120
Flow Line Diagram	122
Appendix B Miscellaneous	123
Cleaning Materials	123
EASY-nLC II Consumable Parts List	123
Replacing the Main Power Fuse	124
Installing and Using the RePlay External Device	125
Declaration of Contamination	133
Transport Instructions	134
Index	135

Preface

This guide describes how to use the EASY-nLC™ II nano-flow liquid chromatography (LC) system with software version 2.8. The information provided also applies to EASY-nLC systems that have been upgraded to version 2.8 (regardless of autosampler version). This guide is intended for laboratory technicians using the EASY-nLC high performance liquid chromatography (HPLC) system to execute analytical runs. It assumes basic knowledge of how to use menu-driven software, and familiarity with standard laboratory practices as well as HPLC terminology and practices.

Before you turn on the instrument, Thermo Fisher Scientific recommends reading [Chapter 1, “Introduction,”](#) [Chapter 2, “Installation,”](#) and [Chapter 3, “Instrument Control Software.”](#) While the instrument is running, review [Chapter 4, “Configuring the EASY-nLC II System,”](#) [Chapter 5, “Preparation for Use,”](#) and [Chapter 6, “Running the First Sample.”](#)

Contents

- [Related Documentation](#)
- [Safety and Special Notices](#)
- [Contacting Us](#)
- [Declaration of Conformity](#)

❖ To suggest changes to documentation or to Help

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



Related Documentation

In addition to this guide, Thermo Fisher Scientific provides the following documents for the EASY-nLC instrument:

- *EASY-nLC Virtual Instrument for Xcalibur Quick Reference*
- *EASY-nLC II Preinstallation Requirements Guide*

Safety and Special Notices

Make sure to follow the safety practices presented in this guide, and observe the safety and special notices that appear in boxes.

Safety Practices

The following safety practices will help to ensure the safe operation of your instrument:

- Before removing the protective panels, disconnect the HPLC from all power sources, or risk exposure to dangerous voltages.
- Always replace blown fuses with fuses of the size and rating indicated on the fuse panel. See [“Replacing the Main Power Fuse.”](#)
- Repair or replace faulty insulation on power cords.
- Use the EASY-nLC II instrument with appliances and power sources that have the proper protective grounding.

When connected to analytical equipment such as a mass spectrometer, also ensure that the protective grounding is shared between the instruments.

- Observe all written safety precautions during all phases of operation, service, and repair of this instrument. Failure to comply with these precautions or with specific warnings elsewhere in this manual violates safety standards of design, manufacture, and intended use of the instrument and may result in injury or loss of life.

Special Notices

The following two stickers appear on the EASY-nLC II HPLC instrument:



This sticker indicates that care should be taken to prevent personal injury or damage to parts of the EASY-nLC II.



This sticker alerts you to consult this manual for instructions on how to operate the HPLC.

The safety and special notices in the documentation include the following:



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 may 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

Fax	(+45) 6557 2301
Web site	www.proxeon.com/productrange/nano_lc
E-mail	easysupport@proxeon.com
Address	Thermo Fisher Scientific Edisonsvej 4 DK-5000 Odense C

Find software updates and utilities to download at mssupport.thermo.com.

❖ To copy manuals from the Internet

Go to mssupport.thermo.com and click **Customer Manuals** in the left margin of the window.

❖ **To contact Customer Service for ordering information¹**

- For North America, see [Table 1](#).
- For Europe, see [Table 2](#).
- For Australia and Asia, see [Table 3 on page xiii](#).

To order consumable and spare parts, go to www.proxeon.com/productrange/nano_lc/spares_accessories.

Table 1. Customer Service information for North America

North America	
United States	
Phone	[1] 800-532-4752
Fax	[1] 877-373-4006
E-mail	us.customer-support.analyze@thermofisher.com
Address	1400 North Point Pkwy #10 West Palm Beach, FL 33407
Canada	
Phone	[1] 800 530 8447
Fax	[1] (905) 890-9161
E-mail	us.customer-support.analyze@thermofisher.com
Address	2845 Argentia Road, Unit 4 Mississauga, Ontario, L5N 8G6

Table 2. Customer Service information for Europe (Sheet 1 of 4)

Europe	
Austria	
Phone	[43] (0) 1 333 50 34-0
Fax	[43] (0) 1 333 50 34-26
E-mail	service.sid.austria@thermofisher.com
Address	Wehlistrasse 27b A-1200 Wien

¹ In the following tables, the country code is enclosed in square brackets []. The city code or area code is enclosed in parenthesis (). For countries other than the U.S., when you are dialing from within the specified country, dial the 0 of the city code. For countries other than Italy, when you are dialing from outside the country, do not dial the 0 of the city code.

Table 2. Customer Service information for Europe (Sheet 2 of 4)**Belgium**

Phone	[32] (0) 2 482 3030
Fax	[32] (0) 2 482 3031
E-mail	service.sid.benelux@thermofisher.com
Address	Clintonpark “Keppekouter” Ninovesteenweg 198 B-9320 ERMEBODEGEM - AALST

Denmark

Phone	[45] (70) 236267
Fax	[45] (70) 236263
E-mail	service.sid.dk@thermofisher.com
Address	Fruebjergvej 3 2100 København Ø

Finland – see Sweden, Norway, and Finland**France****(Also representing French-speaking North Africa, Algeria, Morocco, and Tunisia)**

Phone	[33] (0) 1 60 92 49 50
Fax	[33] (0) 1 60 92 48 99
E-mail	service.sid.lesulis@thermofisher.com
Address	16 Avenue du Québec Silic 765 Z.A. de Courtaboeuf F-91963 Les Ulis Cédex

Germany

Phone	[49] (0) 6103 408 1050
Fax	[49] (0) 6103 408 1213
E-mail	service.dreieich@thermofisher.com
Address	Im Steingrund 4-6 D-63303 Dreieich

Italy

Phone	Numero Verde (800) 823 162
Fax	[39] (02) 95320 225
E-mail	assistenza.tecnica.it@thermofisher.com
Address	Strada Rivoltana I-20090 Rodano (Milano)

Table 2. Customer Service information for Europe (Sheet 3 of 4)**Netherlands**

Phone	[31] (0) 76 579 55 55
Fax	[31] (0) 76 581 09 61
E-mail	service.sid.benelux@thermofisher.com
Address	Takkebijsters 1 NL-4817 BL Breda

Norway – see Sweden, Norway, and Finland**Spain**

Phone	[34] (914) 845 965
Fax	[34] (914) 843 598
E-mail	service.sid.spain@thermofisher.com
Address	C/Valportillo I, n°22 1a Planta Edificio Caoba ES-28108 Alcobendas - Madrid

Sweden, Norway, and Finland

Phone	[46] (0) 8 556 468 20
Fax	[46] (0) 8 556 468 08
E-mail	service.sid.nordic@thermofisher.com
Address	Pyramidbacken 3 S-141 75 Kungens Kurva (Stockholm) Sweden

Switzerland

Phone	[41] (617) 16 77 40
Fax	[41] (617) 16 77 20
E-mail	service.sid.ch@thermofisher.com
Address	Neuhofstrasse 11 4153 Reinach

Table 2. Customer Service information for Europe (Sheet 4 of 4)**United Kingdom**

Phone	[44] (0) 870 241 1034
Fax	[44] (0) 1442 233 667
E-mail	service.sid.hemel@thermofisher.com
Address	Stafford House 1 Boundary Park Boundary Way Hemel Hempstead Hertfordshire HP2 7GE

Table 3. Customer Service information for Australia and Asia (Sheet 1 of 2)**Australia and Asia****Australia**

Phone	[61] 39757 4300
Fax	[61] 9763 1169
E-mail	analyze.au@thermofisher.com
Address	P.O. Box 9092 5 Caribbean Drive Scoresby, VIC 3179

Japan

Phone	[81] (45) 453-9100
Fax	[81] (45) 453-9110
E-mail	analyze.jp@thermofisher.com
Address	C-2F 3-9 Moriya-cho, Kanagawa-ku Yokohama 221-0022

P.R. China

Phone	(free lines) 800 810 5118, 400 650 5118
Fax	[86] 10 88370548
E-mail	analyze.cn@thermofisher.com
Address	7th Floor, 7F Tower West, Younghe Plaza No. 28, Andingmen East Street Beijing 100007

Table 3. Customer Service information for Australia and Asia (Sheet 2 of 2)

P.R. China

Phone	(free lines) 800 810 5118, 400 650 5118
E-mail	analyze.cn@thermofisher.com
Address	Building 6 No. 27 Xin Jin Qiao Road Shanghai 201206

Declaration of Conformity

U.S. Safety and EMC (Electromagnetic Compliance) Standards

Safety

This instrument has been reviewed for compliance with standard ANSI/UL 3101-1, “Electrical Equipment for Laboratory Use; Part 1: General Requirements,” 1st Edition.

EMC

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 undesired operation.



WARNING Changes or modification to this unit not expressly approved by the party responsible for compliance could void the user’s authority to operate the equipment.

Note

This equipment has been tested and found to comply with the limits for a Class A digital device, pursuant to Part 15 of the FCC Rules. These limits are designed to provide reasonable protection against harmful interference when the equipment is operated in a commercial environment. This equipment generates, uses, and can radiate radio frequency energy and, if not installed and used in accordance with the instruction manual, may cause harmful interference to radio communications. Operation of this equipment in a residential area is likely to cause harmful interference in which case the user will be required to correct the interference at his own expense.

Note You must use shielded cables with this unit to ensure compliance with the Class A FCC limits.

Canadian Safety and EMC (Electromagnetic Compliance) Standards

Safety

This instrument has been reviewed for compliance with standard CAN/CSA-C22.2 No. 61010-1, Second edition - "Safety Requirements for Electrical Equipment for Measurement, Control, and Laboratory Use; Part 1: General Requirements."

EMC

This Class A digital apparatus meets all requirements of the Canadian Interference-Causing Equipment Regulations.

European Safety and EMC (Electromagnetic Compliance) Standards

Application of Council Directive(s)	2006/95/EEC "Low Voltage": Intertek Group plc 89/336/EEC "Electromagnetic Compatibility": DELTA Denmark
Standard(s) to which conformity is declared	EN61010-1:2001, Second edition - "Safety Requirements for Electrical Equipment for Measurement, Control and Laboratory Use" EN61010-2-81 "Part 2-081: Particular requirements for automatic and semi-automatic laboratory equipment for analysis and other purposes" EN/(IEC) 61326-1:2006, A1(1998), A2(2001) and A3(2003) "EMC requirements for electrical equipment for measurement, control and laboratory use"
Manufacturer's Name	Proxeon Biosystems A/S
Manufacturer's Address	Edisonsvej 4, DK-5000 Odense, Denmark
Type of Equipment	Laboratory Instrumentation
Model Name	EASY-nLC™ II
Model Numbers	LC110/LC111
Serial Number	LC-000100 and later
Year of Manufacture	2010–

I, the undersigned hereby declare that the equipment specified above conforms to the above Directive(s) and Standard(s).



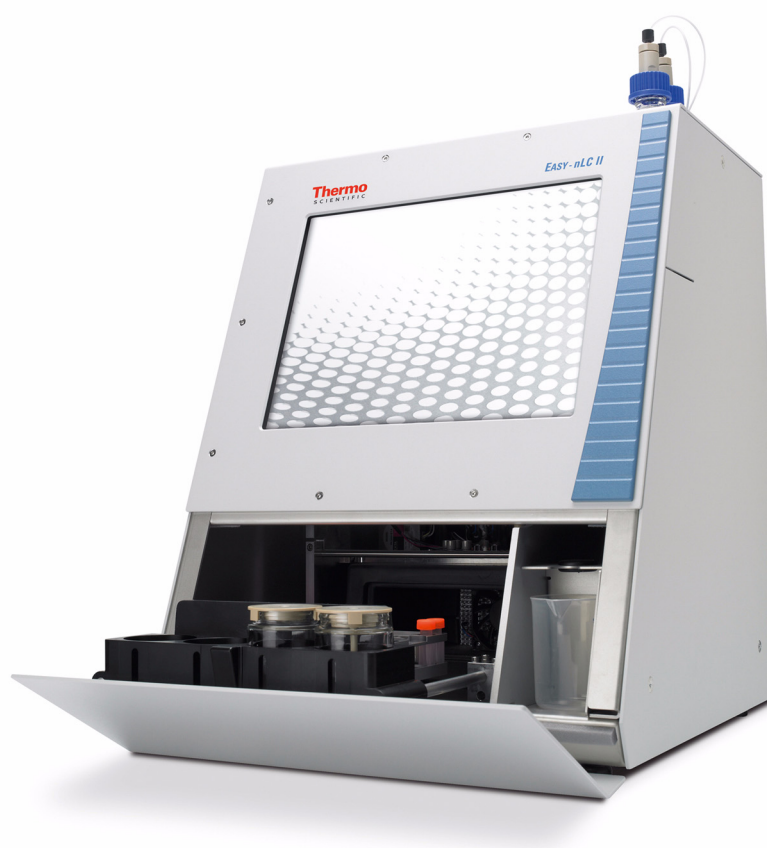
Ole Vorm, Site Manager

August 31, 2010

Introduction

The EASY-nLC II is a compact, high performance liquid chromatography system designed for split-free flows, down to the low nanoLiter/min range. The system provides ease of use and ease of service without sacrificing functionality or performance.

Figure 1. EASY-nLC II system



Contents

- [Product Description](#)
- [Graphical Hardware Overview](#)

Product Description

In the EASY-nLC II system, all necessary components are integrated into one simple, compact unit, making installation easy: connect the power cable and a network cable (if needed) and turn on the power. An embedded computer controls the system and provides an intuitive HPLC interface using an integrated touch-sensitive monitor.

The instrument has a fixed configuration that optimizes the sample handling by using in-depth knowledge of the instrument design and performance characteristics. Instead of having to set up every aspect of a method—from timing to individual control commands—you can enter the parameters that describe the result you want. The software manages the required low-level control.

Furthermore, you can connect the instrument and the integrated computer to the major mass spectrometers and include them in IT schedules for backup procedures and other functions. System maintenance is provided through a range of controlled software, built-in diagnostic tests, and the internal layout of all instrument components.

System Components

The EASY-nLC II system consists of the following components.

Table 1. EASY-nLC II system components (Sheet 1 of 2)

Hardware	
• Pumps	3 high-pressure pumps (1–300 bar) with associated pressure sensors
• Valves	4 high-pressure valves
• Flow sensor	2 nano-flow sensors
• Autosampler	ASC (latest model): <ul style="list-style-type: none">• Cooled (maximum 20 °C [68 °F] below ambient temperature)• Can hold 6 fixed vials, plus either one (1) Microtiter™ plate (96 or 384 wells) or one (1) 6 × 8 vial adapter (for a total of 54 vials)• 4 liquid containers for solvents, rinse, and waste ASA (previous model): <ul style="list-style-type: none">• Cooled (maximum 20 °C [68 °F] below ambient temperature)• Can hold a 4 × 6 vial adapter or one (1) Microtiter plate with 96 or 384 wells• 4 liquid containers for solvents, rinse, and waste
• Computer	Embedded computer

Table 1. EASY-nLC II system components (Sheet 2 of 2)

- Monitor Touch sensitive 12.1 in. (30.7 cm, 800 × 600 pixels) that can be used with gloves

Total size: *w* 35 cm (14 in.), *d* 38 cm (15 in.), *h* 45 cm (18 in.)

Total weight: 32 kg (71 lb)

Utilities: Main power, 120/230 Vac, 50/60 Hz, 250 W

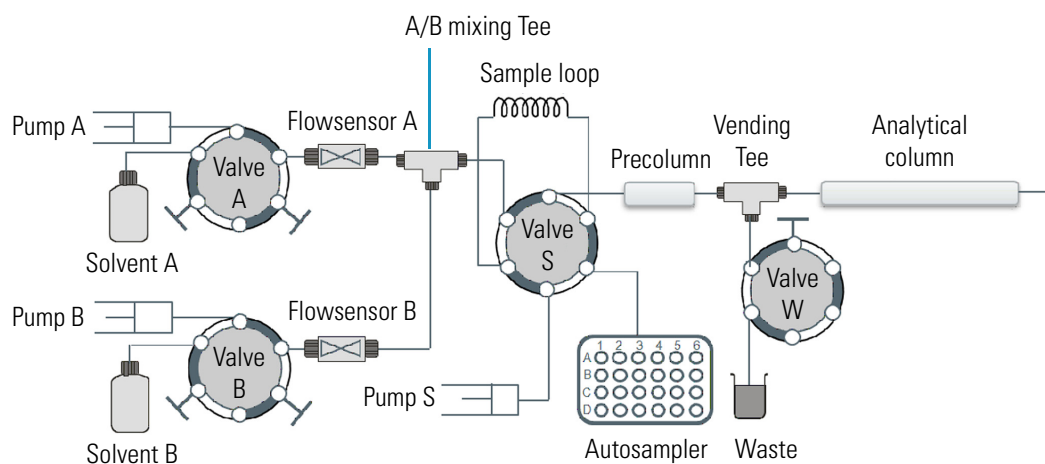
Software

- EASY-nLC II stand-alone application (version 2.8)

System Configuration for Two-Column Setup

For a two-column setup, the components are connected as shown in [Figure 2](#).

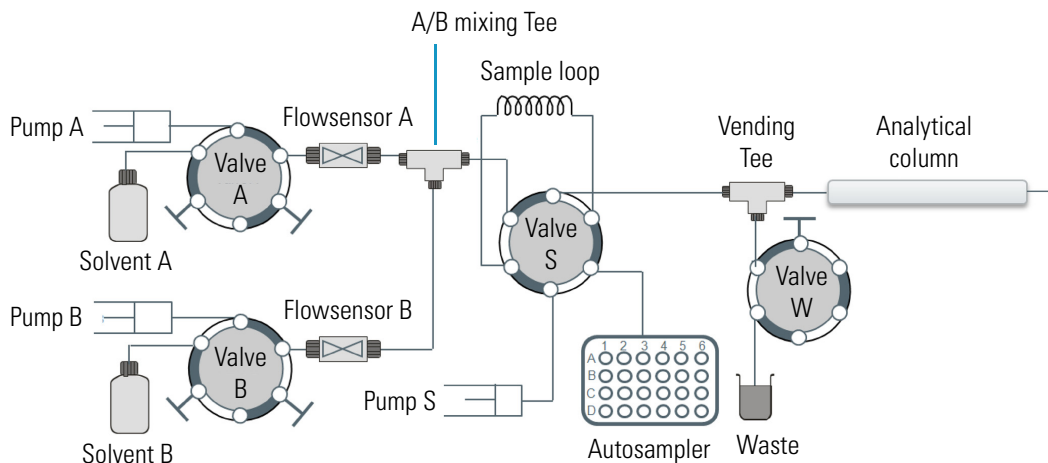
Figure 2. HPLC flow paths for a two-column setup



System Configuration for a One-Column Setup

For a one-column setup, the components are connected as shown in [Figure 3](#).

Figure 3. HPLC flow paths for a one-column setup



Instrument Operation

During sample runs (the execution of methods), the EASY-nLC II instrument goes through a series of predefined steps:

- **Pump refilling** (from solvent bottles)
- **Column reequilibration** (using the A pump)
- **Sample pickup** (where sample is drawn from the autosampler and into the loop)
- **Sample loading** (when the A pump loads the sample from the loop onto the column)
- **Gradient** (formed by the A and B pumps, the sample content is eluted off of the column)
- **Autosampler wash** (when the S pump washes the autosampler needle and loop)

The system performs as many steps in parallel as is physically possible, minimizing the cycle time.

Pump Flow Control

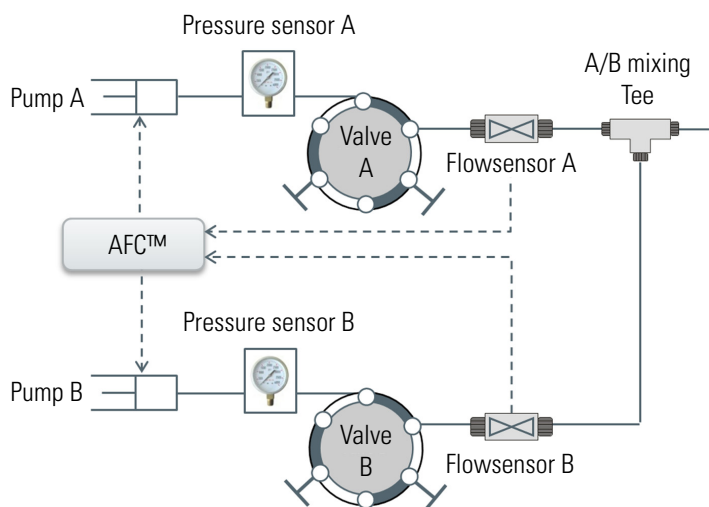
The EASY-nLC II instrument contains split-free, high-pressure syringe pumps capable of delivering flows between 10 nL/min and 300 μ L/min (at different resolutions and with different precision). Different flow control systems are active during different method subprocesses (reequilibration, loading, and gradient) to optimize the pumping system performance. The Advanced Flow Control system (AFC™) accurately controls the flow from each pump (A and B pump), providing enhanced gradient control and retention time reproducibility.

During column reequilibration and sample loading, it is more important to actively control the pressure that these processes are performed under than to have precise flow control (needed during gradient elution). To meet this need, you can use the Intelligent Flow Control system (IFC™) to specify an operating pressure rather than a flow.

Advanced Flow Control System

The advanced flow control system is active during the gradient step of the method execution (and during execution of the isocratic flow maintenance script; see “Prepare – Isocratic Flow” on page 110). The AFC system uses the output of the flow sensors to regulate the flow, accurately maintaining the desired flow from each channel, even during sudden pressure changes that can occur when changing the solvent composition. See Figure 4.

Figure 4. Schematic diagram of the AFC principle

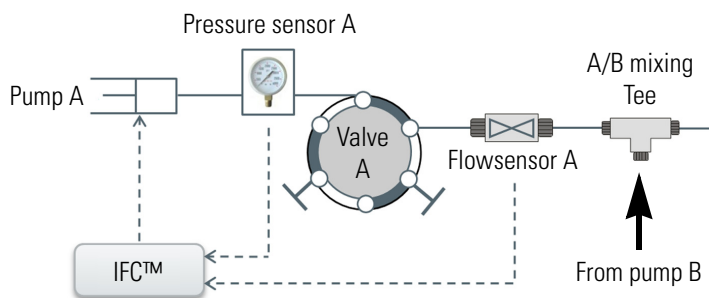


Intelligent Flow Control System

The Intelligent Flow Control (IFC) system is active during reequilibration steps and sample loading, during the execution of the precolumn and analytical column reequilibration scripts (see “Prepare – Precolumn Equilibration” on page 108), and during pressure buildup just before the gradient proper.

The IFC system uses output from both the flow sensor and the pressure sensor to regulate the flow (see Figure 5). You can specify a desired flow rate or a desired operating pressure. This ability reduces the loss of productivity caused by system stoppage related to overpressure (for instance, stoppage caused by bad samples). It also more effectively uses the pressure range of the system (0–300 bar). Additionally, IFC builds up pressure rapidly.

Figure 5. Schematic diagram illustrating the IFC principle



For example, you can program the system to reequilibrate columns at a pressure of 180 bar, at 5 $\mu\text{L}/\text{min}$, or at both of these settings. The setting reached first (which becomes the limiting parameter) depends on the dimensions of the attached column or columns. If the instrument reaches 180 bar pressure before it reaches the 5 $\mu\text{L}/\text{min}$ set point, it adjusts the flow rate to maintain the 180 bar pressure. If the 5 $\mu\text{L}/\text{min}$ is reached before the 180 bar pressure, the pump adjusts the flow rate to maintain the 5 $\mu\text{L}/\text{min}$ through the flow sensor. The maximum flow with the IFC algorithm is 100 $\mu\text{L}/\text{min}$, but the back pressure of the system itself, without columns attached, reaches maximum pressure between 20–25 $\mu\text{L}/\text{min}$.

The duration of the reequilibration or loading step depends on the total volume of solvent specified and the actual flow rate. If the set pressure is the limiting factor, the flow rate is variable as a consequence, varying the duration of these steps. If the duration of the reequilibration and loading steps must be fixed, limit the step by flow rather than the pressure.

Using the EASY-nLC II application, you can turn off the IFC system for reequilibration and loading. The EASY-nLC system then performs these two steps, ignoring any input in the Max Pressure input box of the method editor and regulating the pumps passively by using the specified flow rate.

To reduce cycle time and aid system robustness, keep the IFC system turned on.

Graphical Hardware Overview

The following figures illustrate both external and internal features of the EASY-nLC II instrument. [Table 2](#) describes the elements on the instrument's back panel.

Figure 6. External system features

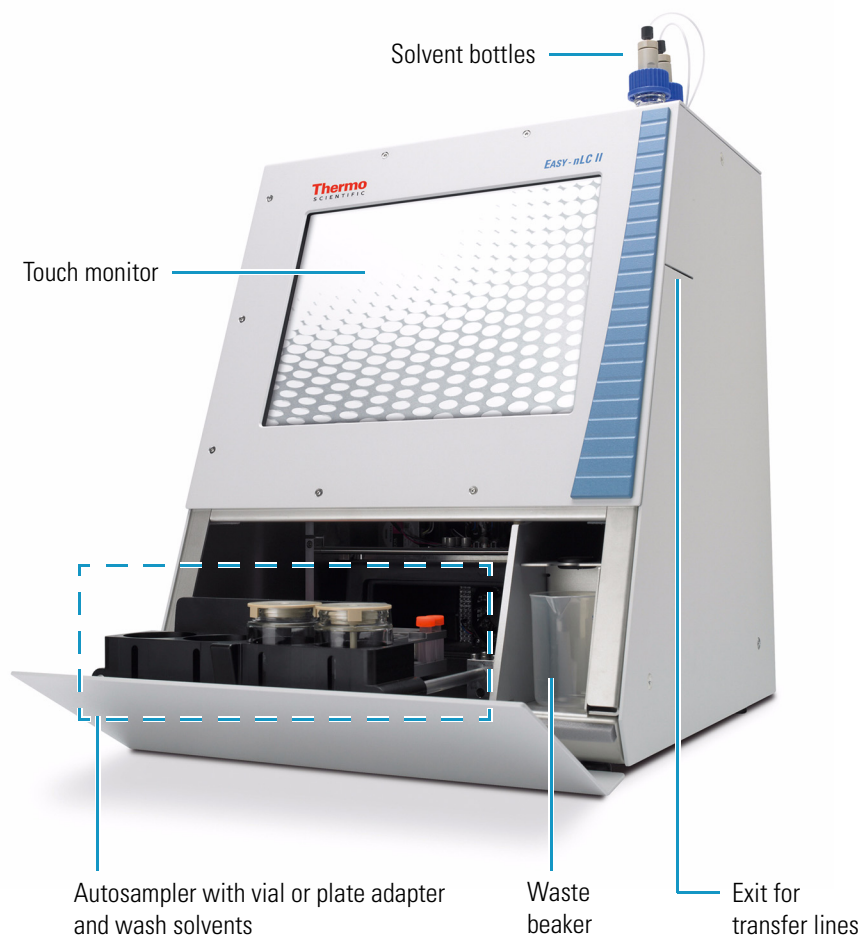


Figure 7. External view of instrument's back panel

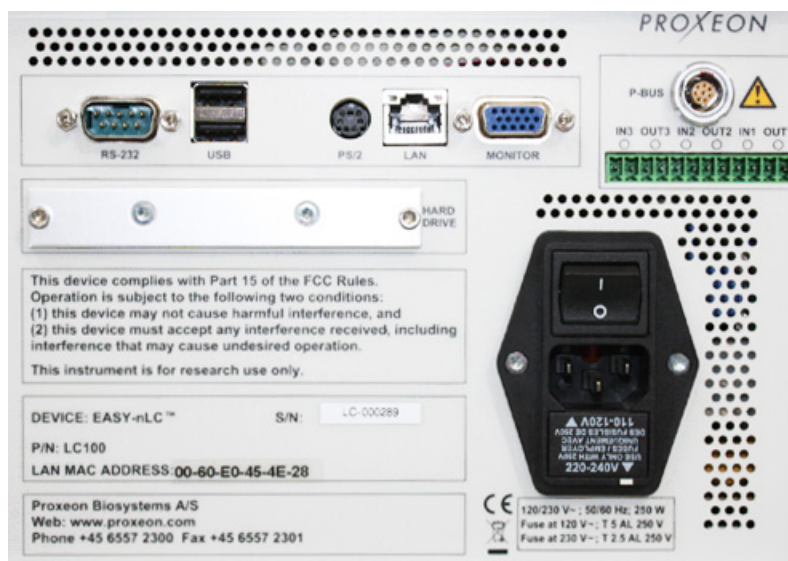


Table 2. Back panel elements

Element	Description
IN/OUT	Contact closure (primarily for communication with the mass spectrometer)
P-BUS	For communication with HPLC add-on instruments
RS-232	For communication with HPLC add-on instruments, for example a syringe drive
USB	Port for keyboard or mouse, two (2) provided
PS/2	Input for connection of keyboard and mouse
LAN	For 10/100 Mb/sec Ethernet connection
MONITOR	Output for connection of external display
I O	Mains power switch
Fuse holder below I O switch	Replaceable fuse ratings: <ul style="list-style-type: none"> • For 120 V, T 5 AL, 250 V • For 230 V, T 2.5 AL, 250 V
P/N	Product order code
S/N	HPLC serial number
LAN MAC ADDRESS	The MAC address of the embedded computer (Gives the EASY-nLC II system a fixed IP address on your local network.)

Figure 8. Internal features (view 1)

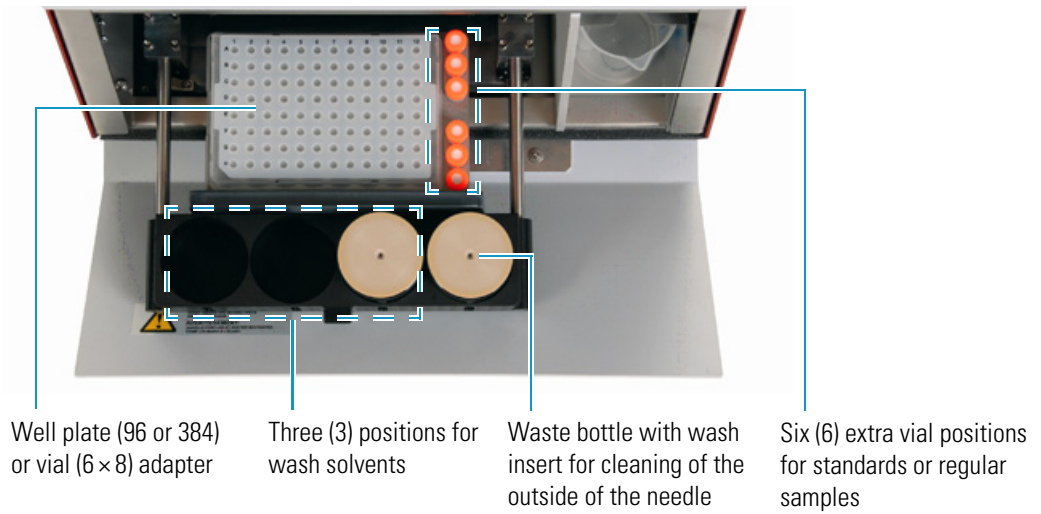


Figure 9. Internal features (view 2)

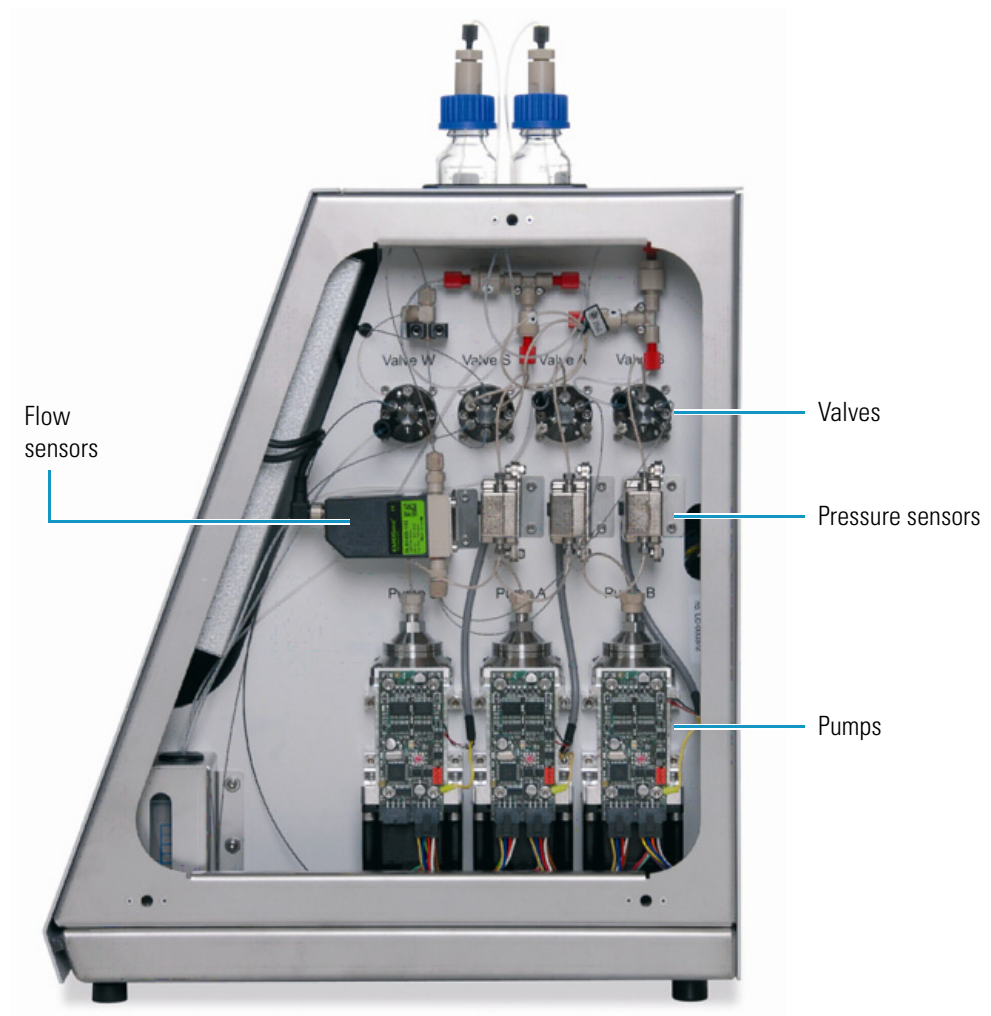
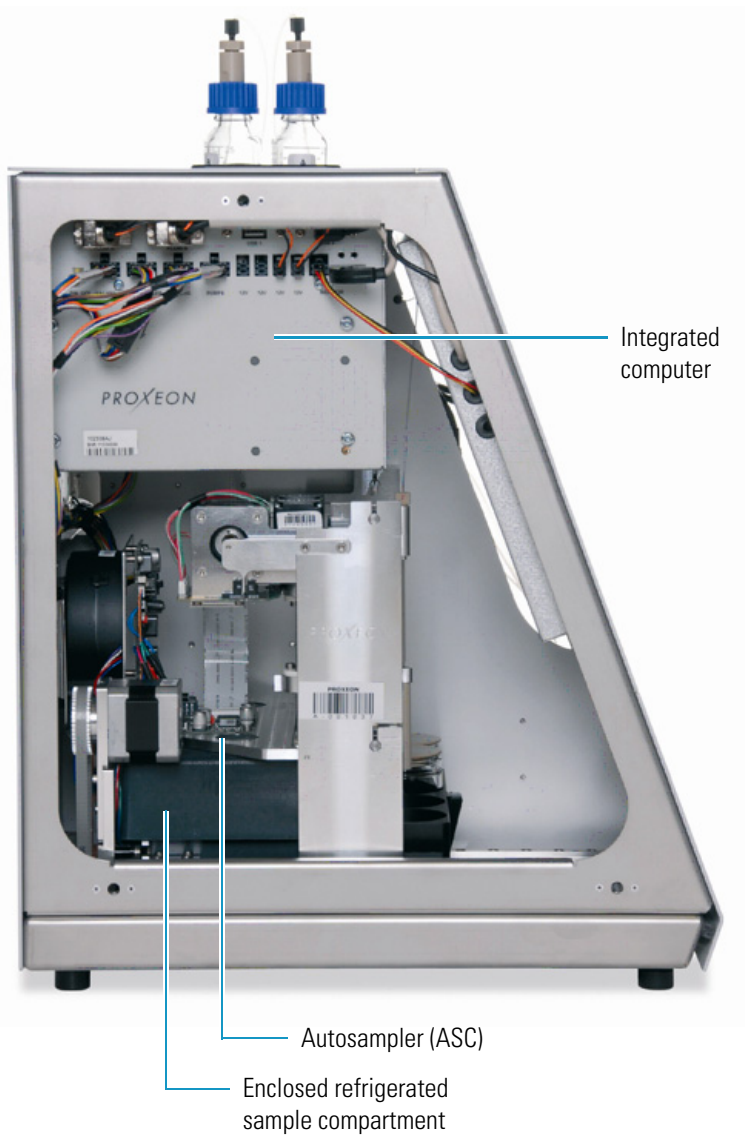


Figure 10. Internal features (view 3)



Installation

The following sections provide important information on installing the Thermo Scientific EASY-nLC II instrument. Be sure to review the “[Safety and Special Notices](#)” on page viii before proceeding.

Contents

- [Placement](#)
- [Power](#)
- [Temperature and Humidity](#)
- [Lifting Instruction](#)
- [External Connections](#)
- [Flow Parts and Lines](#)

Placement

The EASY-nLC II instrument weighs approximately 32 kg (71 lb) and its dimensions are approximately $w 35 \times d 38 \times h 45$ cm ($w 14 \times d 15 \times h 18$ in.). Allow for at least 15 cm (6 in.) of free space at the back of the instrument for proper air circulation.

Power

Use this instrument only with properly grounded appliances and power sources.

120 Vac, 50/60 Hz, 250 W

230 Vac, 50/60 Hz, 250 W

(For UPS dimensioning, assume 250 W)

For 120 Vac: one T 5 AL 250 V fuse (5 × 20 mm, IEC 60127)

For 230 Vac: one T 2.5 AL 250 V fuse

All fuses supplied with the instrument are UL Listed and CSA-certified.

Temperature and Humidity

Avoid locations with high air humidity or extreme changes in temperature (such as direct sunlight, drafts, or directly below an air conditioning or heating vent). Place the EASY-nLC II instrument as close as possible to the detector and mass spectrometer to minimize dead volume in the transfer lines. Place the EASY-nLC II system away from any detector/mass spectrometer vents, however, to avoid exposure to direct heat.

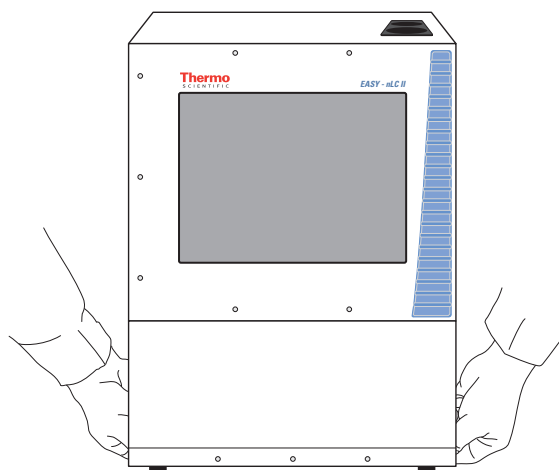
For optimal autosampler plate cooling, place the unit in an area where the working temperature is within 5–30 °C (41–86 °F) .

The optimal humidity range is between 20–80% RH. Avoid condensing humidity.

Lifting Instruction

The instrument weights 32 kg (71 lb). For safety reasons, use two people to move the instrument to a table cart for transport. Gloves are recommended.

Figure 11. Lifting by hand



External Connections

The EASY-nLC II instrument has several types of connections to external services or devices that you can make:

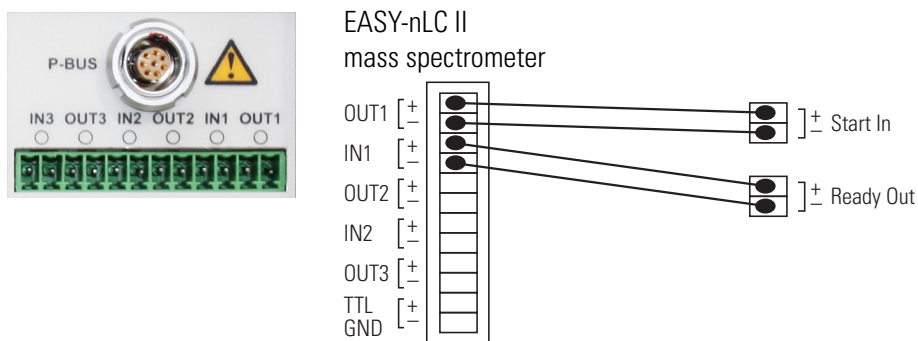
- [Connecting to the Mass Spectrometer through Contact Closure](#)
- [Connecting to Remote Support](#)
- [Connecting to a Local Area Network for Data Exchange](#)
- [Attaching Mouse and Keyboard to the USB Connections](#)
- [Attaching Add-on Products through the P-Bus and RS-232](#)

Connecting to the Mass Spectrometer through Contact Closure

The EASY-nLC II stand-alone application contains preconfigured values for most of the standard mass spectrometers used in proteomics. For instructions on setting up the mass spectrometer connection in the application, see “[Setting Up the Mass Spectrometer Connection](#)” on [page 35](#).

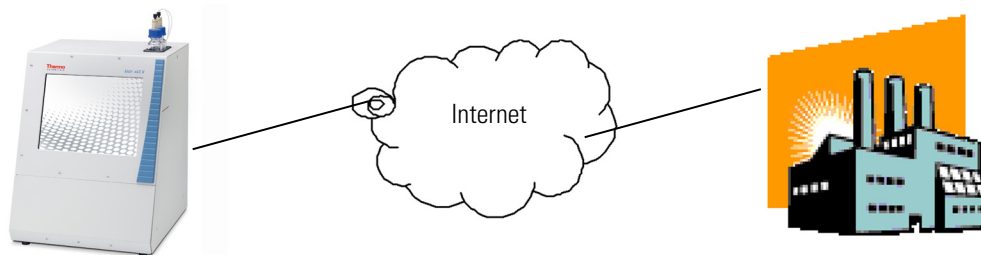
To connect a mass spectrometer, see [Figure 12](#).

Figure 12. Two-way plug configuration for a mass spectrometer



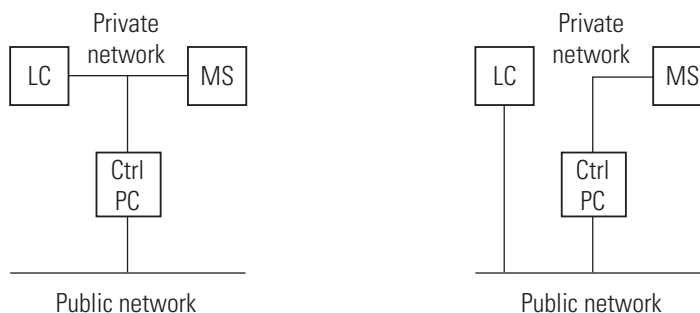
Connecting to Remote Support

You can have Thermo Fisher Scientific online support review your EASY-nLC II system. Using shielded cables, connect the LAN port of the instrument to the local network that has Internet capability.



If the EASY-nLC II system sits on the private network side of the mass spectrometer (MS) control computer, you must move it to the public network, permanently or temporarily, to access the remote support server. See [Figure 13](#).

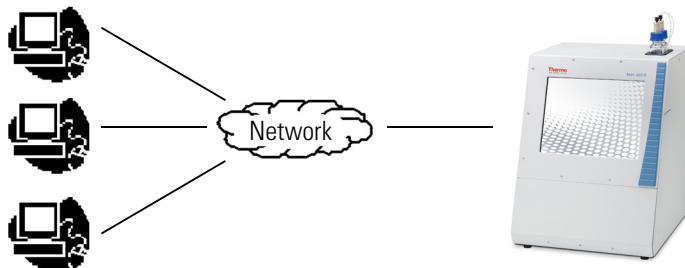
Figure 13. Mass spectrometer control computer from private to public network



Connecting to a Local Area Network for Data Exchange

The EASY-nLC II instrument can work in stand-alone mode, but if you require access to a network, ask the local network administrator for guidance (see “[Network Connection](#)” on page 36).

Figure 14. Connecting the HPLC to a computer network



The instrument can be connected to ordinary computer networks by using a standard Ethernet cable that has RJ-45 connectors. When the instrument connects to a network, other computers on the network can access files on the instrument (for example, logs, exports, imports, reports, statistics, and backups) when they know the IP address of the instrument and have a valid user login account and password.

Attaching Mouse and Keyboard to the USB Connections

You can use a mouse and keyboard instead of the built-in touch screen, but Thermo Fisher Scientific recommends using the touch interface. For more information, see “[Using the Touch-Sensitive Screen](#)” on page 18.

Connect the USB mouse or keyboard by using the USB connectors on the back of the instrument. The EASY-nLC II application detects the mouse or keyboard within 30 seconds. The application does not recognize all USB-based keyboards or mice.

Attaching Add-on Products through the P-Bus and RS-232

The add-ons are controlled and powered through the P-bus and/or controlled through the RS-232 interface.

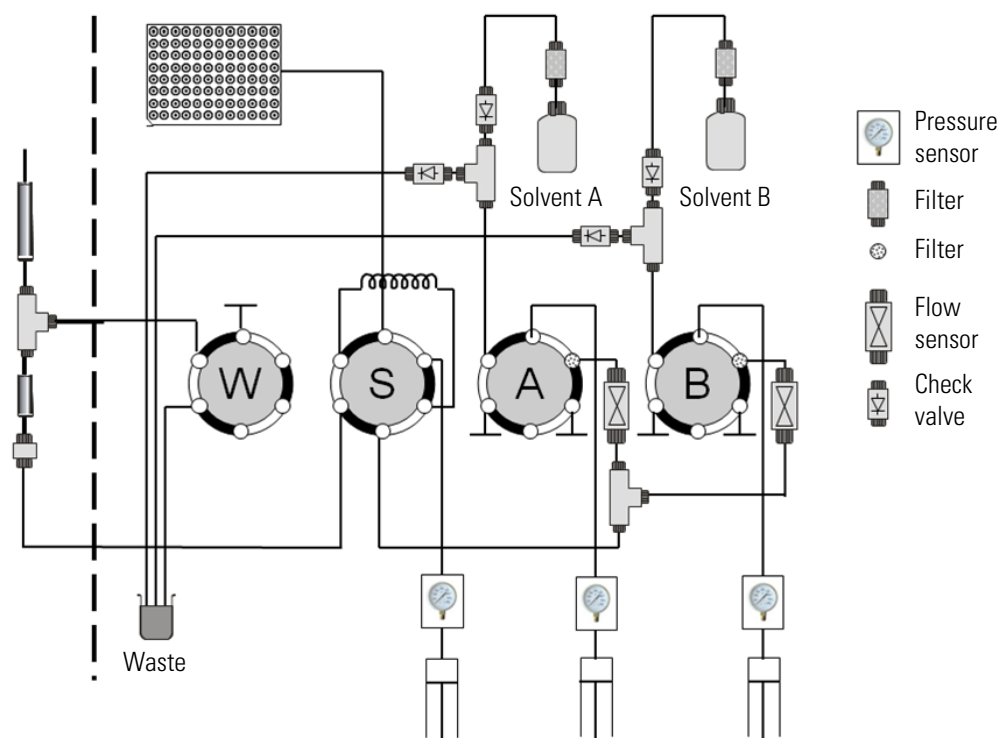


Flow Parts and Lines

The pump and valve compartment of the EASY-nLC II instrument also contains all the flow lines between the components (see [Figure 15](#)).

In addition to pumps and valves, several filters and check valves have been added. The two check valves on lines A and B ensure that solvent is drawn from the solvent bottles and ejected to waste without switching valve positions. The filters and check valves are not visible on most schematics because these extra components do not affect the main sample processing logic.

Figure 15. Complete physical layout and schematics



For complete dimensions, see [“Technical Specifications”](#) on [page 120](#).

Instrument Control Software

The EASY-nLC II instrument is controlled by software that runs on an integrated computer; this eliminates the need for a dedicated computer next to the instrument. This chapter provides basic instructions for the reliable and efficient operation of the EASY-nLC II system.

The application presents an easy-to-use interface for defining methods (sample processing protocols), scheduling batches for execution (on a Microtiter-plate basis), following the system progress and status, managing users, and carrying out maintenance and repair work.

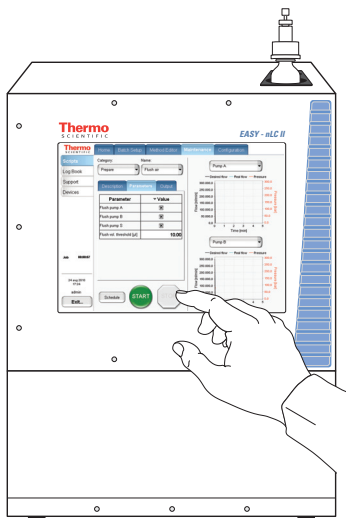
Contents

- [Using the Touch-Sensitive Screen](#)
- [Starting the EASY-nLC II System](#)
- [User Interface Layout](#)
- [General Interaction Principles](#)
- [Application Menu Structure](#)
- [Logging In to the EASY-nLC II System](#)
- [Closing Down the EASY-nLC II System](#)

Using the Touch-Sensitive Screen

The EASY-nLC II system has no keyboard or computer mouse. Control of the instrument comes directly through the instrument monitor.

Figure 16. Finger-touch control



The monitor face is a transparent layer of glass that detects finger pressure (even while wearing gloves) and sends the corresponding commands to the computer.

In most cases, pressing a button displayed on the screen causes the instrument to carry out a certain operation. However, when the application requires text or numeric input, the application displays a keyboard or numeric keypad on top of the main window (Figure 17) so that you can enter values on the screen by pressing the appropriate keys.

Figure 17. Touch keyboards



The keys light up when pressed so that you do not need to look at the text or number field where the entry is also shown.

If preferred, you can attach a USB-based mouse or keyboard to the instrument's back panel, but the user interface has been optimized for touch-screen interaction.

Starting the EASY-nLC II System

Press the On/Off switch on the back of the instrument.

Figure 18. The On/Off switch on the back panel



The start-up procedure takes between 5 and 10 minutes while the instrument does the following:

1. Boots the internal computer.
2. Initializes all of the necessary software components.
3. Checks all the hardware parts (pumps, valves, autosampler, plate-cooler, flow sensor, and so on).
4. Starts the application.

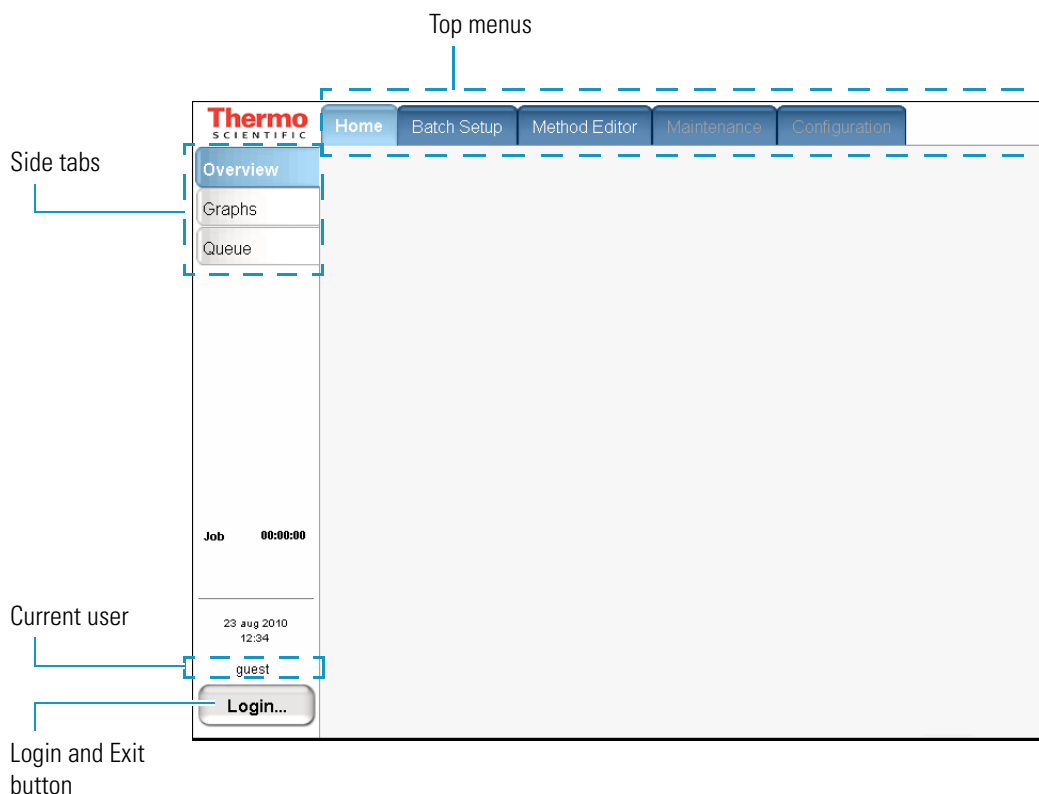
During this process, the screen shows a status bar to indicate progress.



User Interface Layout

The EASY-nLC II user interface provides the following menu structure:

Figure 19. EASY-nLC II user interface layout with menus



Begin by choosing one of the top menus. The application displays a number of tabs down the left side of the active area. Press these tabs to display views related to that topic. During a session, the application remembers which side tab you selected inside each top menu and shows that tab when you next choose that menu.

Use the current user area to log off any page in the application.

Press the logo in the upper left corner to view version information about the currently running software.

General Interaction Principles

Press buttons only one time for a specific action. If the action is important or irreversible, the EASY-nLC II application displays a confirmation dialog box so that you can cancel the action.

You can edit many tables (usually when you have entered the data yourself). You can also access single cells by touching the cell twice to open a keyboard/keypad display for changing the entry.

Long tables have vertical scroll bars that you move either by pressing the up or down arrow symbols, or by dragging the scroll bar and moving it explicitly.

In some tables you can select multiple rows by pressing check boxes on several rows or complete columns by pressing the table headings.

If a screen has blank input fields, press inside the field to enter data. The application displays a keyboard/keypad.

Many button actions provide a confirmation box where the user is prompted to either accept or cancel an action.

In the rest of this guide, we use the menu/left side area to refer to a given screen, accessed first by pressing the top menu tab and then by pressing the appropriate side tab (for example **Home > Queue**).

Application Menu Structure

The EASY-nLC II application has five top menus:

- [Home](#)
- [Batch Setup](#)
- [Method Editor](#)
- [Maintenance](#)
- [Configuration](#)

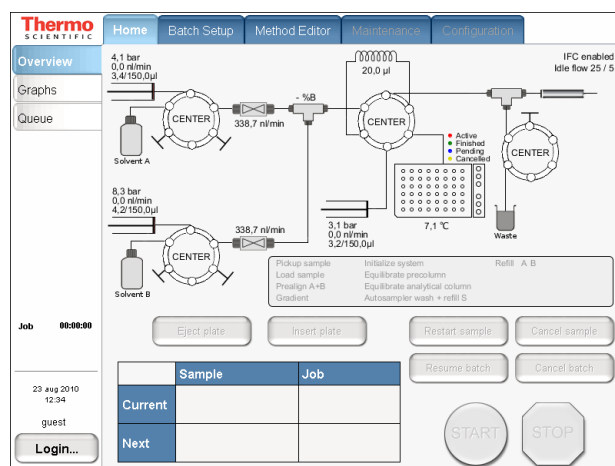
Home

Press the **Home** tab to view the default menu for the application. It provides three pages to view and change information about running samples and instrument components.

Table 3. Home menu pages (Sheet 1 of 2)

Use the **Overview** page to view a graphical layout of the instrument components. Graphics of most hardware parts are made touch sensitive, so you can control them through the shortcut menu that appears when you touch a specific part. This view is dynamically updated.

For more information, see “Using the Overview Page in the Home Menu” on page 64.



Use the **Graphs** page to view up to four graphs (each selectable from a list) with historic data for the currently active sample.

For more information, see “Using the Graphs Page in the Home Menu” on page 65.

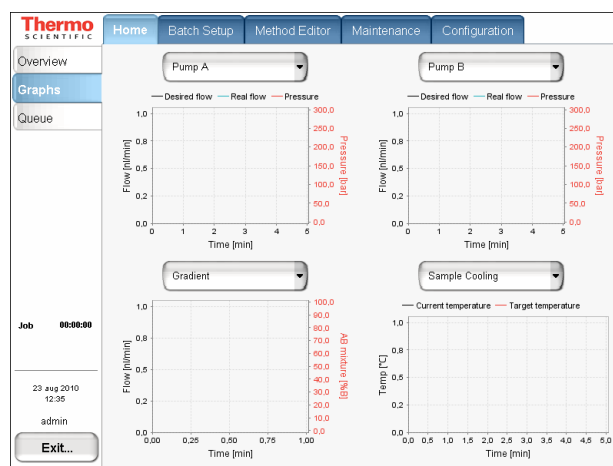
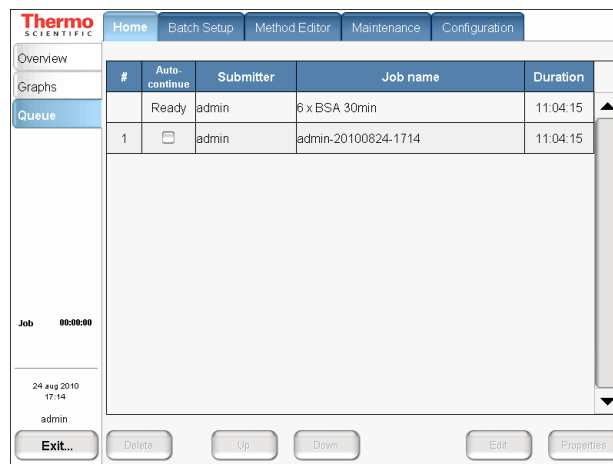


Table 3. Home menu pages (Sheet 2 of 2)

Use the **Queue** page to view a list of batch jobs, currently executing or waiting. You can edit a running batch from this page.

For more information, see “[Editing the Running Batch](#)” on page 67.



Batch Setup

Specify and schedule sample processing jobs for execution by using views under the Batch Setup tab.

Table 4. Batch Setup pages (Sheet 1 of 2)

Use the **File** page to view or change a directory tree of saved batch jobs that can be copied, deleted, or exported. You can create or import new jobs and create or delete directories on this page.

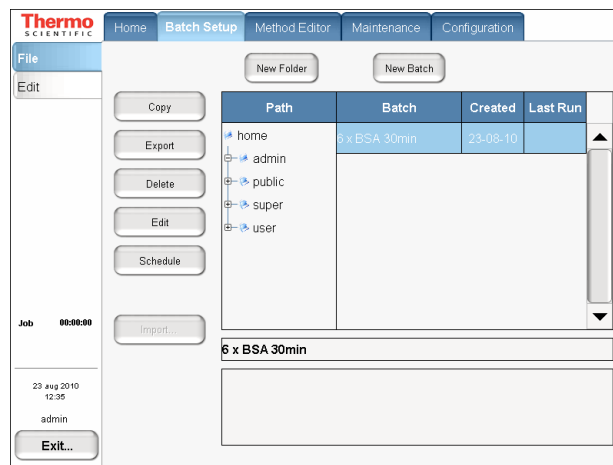
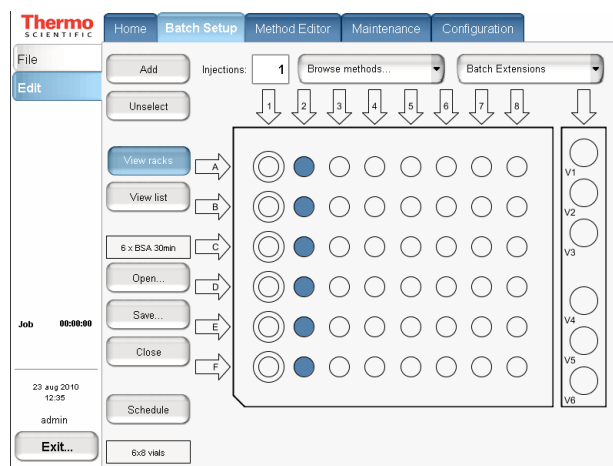


Table 4. Batch Setup pages (Sheet 2 of 2)

Use the **Edit** page to specify new batch jobs. There are two views:

- Use the **View Racks** view to see a graphical view of the chosen plate format. Select a method from the list, enter sample volume and number of injections, highlight the appropriate positions on the plate (use the arrows to select whole rows or columns) and add the selection to the batch job. Repeat the process and either save the batch or schedule it directly.



When using 384 well plates, you can use the arrows near the bottom of the screen to navigate between the currently shown selections. In this case, pressing the middle square shows a full-screen view (not active for input) and pressing anywhere in the full-screen view takes you back to the 8 × 12 active selection.

- Use the **View List** view to see the batch job as a list of samples in the order of processing. Each row in the table holds information about the position, name, and method for that position.

For more information, see [“Editing the Running Batch”](#) on page 67.



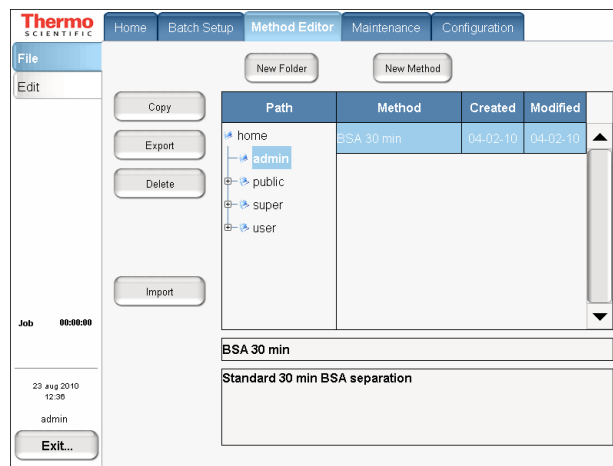
Press **Schedule** to enter the samples for processing. This takes you to the Batch Queue tab in the Home menu.

Method Editor

Press the **Method Editor** tab to define sample processing methods.

Table 5. Method Editor pages (Sheet 1 of 3)

Use the **File** page to view a directory tree of saved methods that you can copy, delete, or export. You can create or import new methods and create or delete directories from this page.



Use the **Edit (1/5)** page to view and change the values that make up a method. The methods are result oriented, so you state the type of performance you require from the instrument. Each section shows the flow parts that are affected or controlled by the inputs.

For more information on this page of the Edit view, see [“Page 1/5”](#) on [page 54](#).

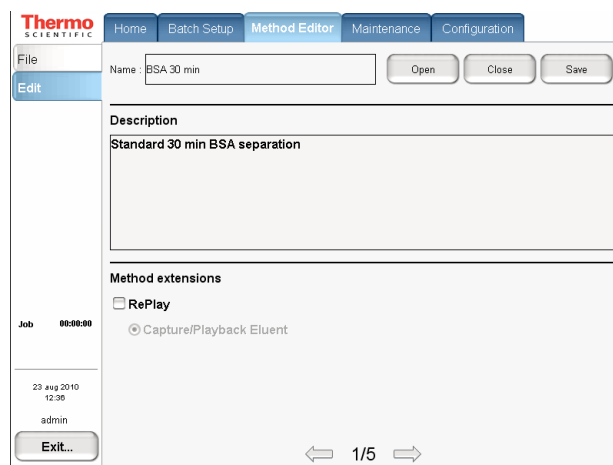


Table 5. Method Editor pages (Sheet 2 of 3)

Use the **Edit (2/5)** page to control sample pickup.

For more information on this page of the Edit view, see “Page 2/5” on page 55.

Use the **Edit (3/5)** page to control the gradient.

For more information on this page of the Edit view, see “Page 3/5” on page 56.

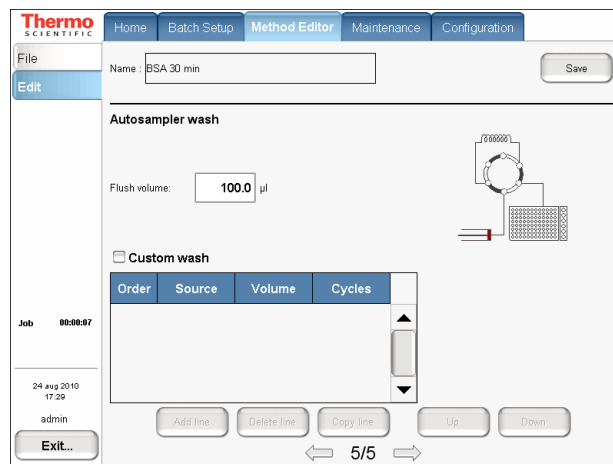
Use the **Edit (4/5)** page to control reequilibration.

For more information on this page of the Edit view, see “Page 4/5” on page 58.

Table 5. Method Editor pages (Sheet 3 of 3)

Use the **Edit (5/5)** page to control autosampler wash.

For more information on this page of the Edit view, see “Page 5/5” on page 59.



Maintenance

Press the **Maintenance** tab to perform a number of functions designed to provide and improve system status.

Table 6. Maintenance pages (Sheet 1 of 2)

Use the **Scripts** page to view a collection of useful maintenance procedures for keeping the instrument functioning properly and help pinpoint possible problems. Select the category and then the individual script. After selection, the application provides explanatory text when the list of input parameters appears. You must enter parameters correctly before starting the script. Next to the script control area, the application displays two configurable graphs so that you can monitor the process.

For the complete list of maintenance scripts and how to set them up, see “Maintenance Scripts” on page 105.

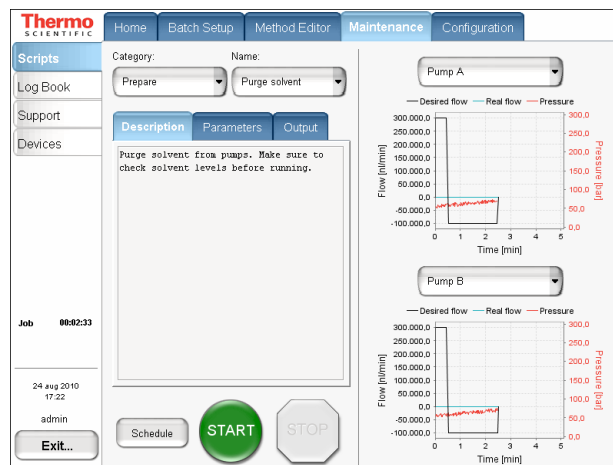
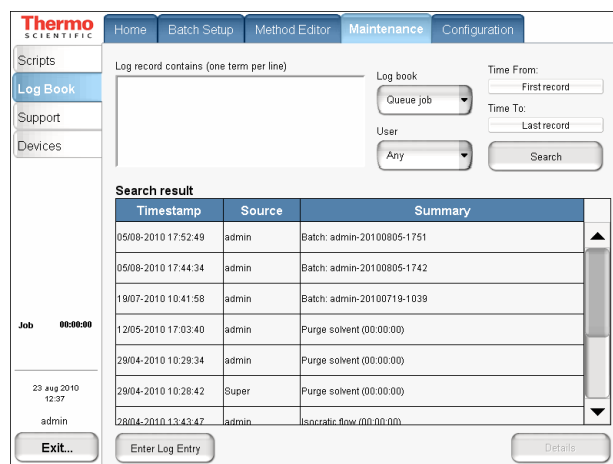
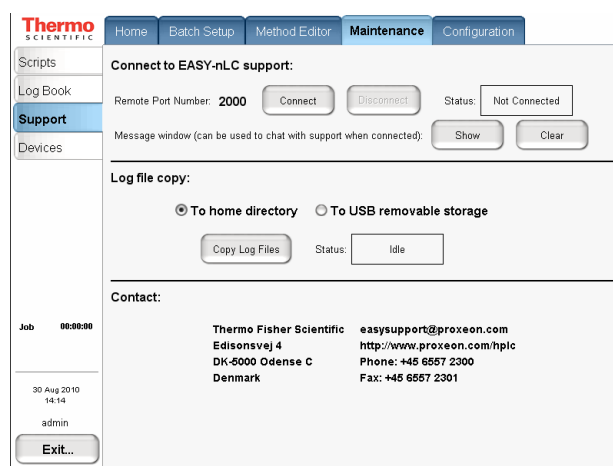


Table 6. Maintenance pages (Sheet 2 of 2)

Use the **Log Book** page to electronically enter all the service actions that have been carried out on the instrument. There is a general log where you can enter general comments about all the components and areas for several specific actions that are used actively in other parts of the application (primarily statistics).

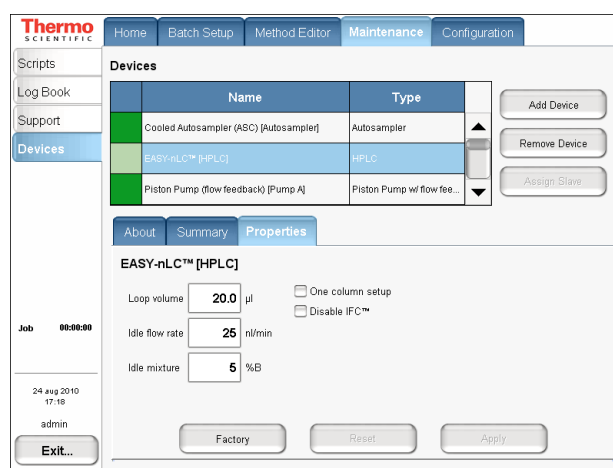


Use the **Support** page to set up support from Thermo Fisher Scientific. For a more comprehensive description, see [Chapter 9, “Remote Support.”](#)



Use the **Devices** page to view and configure hardware devices (including calibrating the autosampler). You can add instruments to the HPLC, such as the Advion™ RePlay™ device (see [“Installing and Using the RePlay External Device”](#) on page 125).

For configuration details and parameter descriptions, see [“Configuring Column Setup, Loop Volume, Idle Flow, and IFC”](#) on page 40



Configuration

Press the **Configuration** tab to control the instrument setup and manage system configuration.

Table 7. Configuration pages (Sheet 1 of 3)

Use the **Users** page to manage user accounts (create, edit, and delete).

For more information, see [“Creating New User Accounts”](#) on page 40.

The screenshot shows the 'Users' configuration page in the Thermo Scientific software. The 'Configuration' tab is selected. On the left, a sidebar contains 'Users', 'Connections', 'Network', 'Time', and 'Backup'. The main area is titled 'Users' and shows a form for editing the 'admin' user. Fields include 'Name' (Administrator), 'Password' (masked with asterisks), and 'Re enter' (masked). A 'Description' field contains 'Administrator'. The 'Level' is set to 'Admin' with radio buttons for 'Normal', 'Super', and 'Admin'. Below are dropdown menus for 'Pump A', 'Pump B', 'Gradient', and 'Sample Cooling'. At the bottom are buttons for 'Open', 'Close', 'Save', 'Delete', and 'New'. A status bar at the bottom left shows 'Job 00:00:00', '24 Aug 2010 17:17', and 'admin'.

Use the **Connections** page to view which mass spectrometer connects to the EASY-nLC II system and how the communication is set up.

For more information, see [“Setting Up the Mass Spectrometer Connection”](#) on page 35.

The screenshot shows the 'Mass Spectrometer' configuration page. The 'Configuration' tab is selected. The sidebar on the left is the same as in the previous screenshot. The main area is titled 'Mass Spectrometer' and shows the 'Instrument' set to 'Thermo Finnigan/LTQ Orbitrap'. Under 'Communications', 'Use Contact Closure' is checked, while 'Use Ethernet' and 'Use RS-232' are unchecked. The 'Contact Closure' section shows settings for 'EASY-nLC™' and 'Thermo Finnigan/LTQ Orbitrap'. The 'Protocol' is 'Two-way', 'State at start' is 'Open', and 'Start at' is 'Gradient'. The 'Signal width' is set to '1 sec.'. A diagram shows connections between the EASY-nLC™ ports (OUT1, IN1, OUT2, IN2, OUT3, TTL, GND) and the Peripheral Control ports (Start In, Ready Out, Ready In, Start Out). Buttons for 'Default', 'Undo Changes', and 'Save' are at the bottom. The status bar at the bottom left shows 'Job 00:00:00', '23 Aug 2010 12:38', and 'admin'.

Table 7. Configuration pages (Sheet 2 of 3)

Use the **Network** page to understand how the EASY-nLC II system is identified in a connected computer network. Consult your IT administrator before making changes to the system.

For more information, see “[Network Connection](#)” on page 36.

The screenshot shows the 'Network' configuration page in the Thermo Scientific software. The interface includes a top navigation bar with 'Home', 'Batch Setup', 'Method Editor', 'Maintenance', and 'Configuration'. A left sidebar contains 'Users', 'Connections', 'Network', 'Time', and 'Backup', with 'Network' selected. The main area is titled 'Address' and features radio buttons for 'Dynamic' (selected) and 'Fixed' network types. Below this, there are input fields for 'Name' (LC-000125) and 'Domain' (proxeon.com). IP address fields are provided for 'IP', 'Subnet', 'Gateway', and 'DNS', each with four digit boxes. A 'Support server' section contains an IP field with values 195, 41, 108, and 93, and a 'Default IP' button. A 'Save configuration' button is located at the bottom right of the main area. The bottom left corner shows 'Job 00:01:02', the date '24 aug 2010 23:42', the user 'admin', and an 'Exit...' button.

Use the **Time** page to set the internal clock in the EASY-nLC II system and the display format used throughout the application to print date and time information.

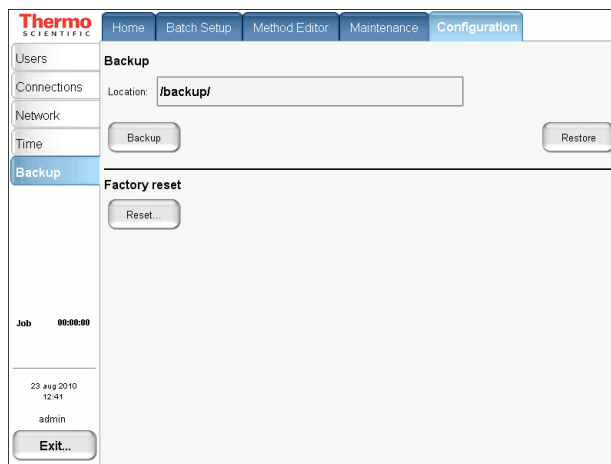
The screenshot shows the 'Time Settings' page in the Thermo Scientific software. The top navigation bar and left sidebar are identical to the previous screenshot, with 'Time' selected. The main area is titled 'Time Settings' and contains several controls: a 'Set Time Zone' dropdown menu with a 'Reset' button; 'Set Time...' and 'Synchronize' buttons; an 'NTP server' field containing 'clock.isc.org'; and a 'Set Format' field showing 'dd/MM-yyyy HH:mm:ss' with a 'Presets' dropdown menu. The bottom left corner displays 'Job 00:00:00', the date '23 aug 2010 12:41', the user 'admin', and an 'Exit...' button.

Table 7. Configuration pages (Sheet 3 of 3)

Use the **Backup** page to save and restore backups of the system database and configuration. A backup includes methods, batches, instrument configuration, and user accounts. Data that you restore from a backup includes all of this information.

You can also remove changes you have made to the system, completely or selectively, by restoring the factory default settings.

Performing a system restoration does not impact statistics on pumps, valves, and so on in the Maintenance/Statistics menu.



IMPORTANT Make sure that no one else is using the EASY-nLC II system when restoring a backup because this also influences user accounts.

Logging In to the EASY-nLC II System

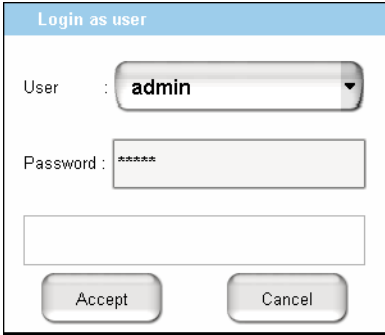
The EASY-nLC II application has a user management system that assigns specific access rights to each user. When the instrument ships, it contains two predefined users: guest and admin. The default user account is “guest” (and does not require a password). With this account you can see and monitor the system, but you cannot enter any new information.

❖ To start using or configuring the EASY-nLC II system

1. Log in as the **admin** user (see [Figure 20](#)).
2. Press **Login** (the factory password for the admin account is also **admin**).

The first time you start the application, it automatically logs you in as a guest. “Guest” appears above the Login button.

Figure 20. EASY-nLC II Login screen

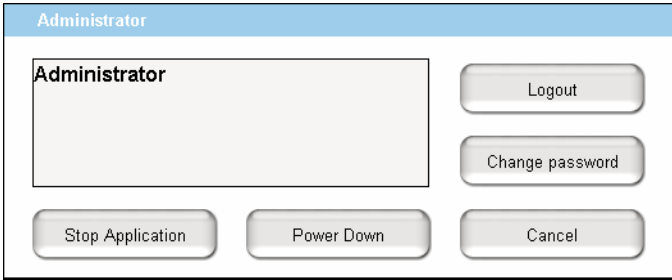


The screenshot shows a dialog box titled "Login as user". It contains a "User" dropdown menu with "admin" selected, a "Password" field with "*****" entered, and an empty field below it. At the bottom are "Accept" and "Cancel" buttons.

IMPORTANT Change the admin password as soon as possible for security reasons.

❖ To log off the EASY-nLC II system

1. Press **Exit**.
2. Press **Logout** in the confirmation dialog box.



The screenshot shows a dialog box titled "Administrator". It contains a large empty box with "Administrator" text, and buttons for "Logout", "Change password", "Stop Application", "Power Down", and "Cancel".

An administrator can start a controlled power-down or exit the application from this dialog box. If you are not an administrator, you can only log out or change your password.

Closing Down the EASY-nLC II System

Closing down the instrument in a controlled manner is important to allow all the components to shut down in an orderly sequence. Using this controlled method saves important data so that the instrument starts with the correct information the next time you use it.

WARNING If you turn off the power switch during normal operation, you risk damaging essential system components. Follow the procedure outlined here whenever possible.

❖ To turn off the EASY-nLC II system

1. Press **Exit**.
2. Press **Power Down**.

The EASY-nLC II application displays a white screen with a small progress bar. When the progress bar is filled out and the message appears indicating you can safely turn off the instrument, go to the next step.

3. After receiving the message that you can safely turn off the instrument, turn off the power switch on the back of the instrument.



Configuring the EASY-nLC II System

After turning the instrument on and logging on as a system administrator, you must configure the instrument for use and define the work environment between the instrument and computer.

Contents

- [Setting Up the Mass Spectrometer Connection](#)
- [Network Connection](#)
- [User Permissions](#)

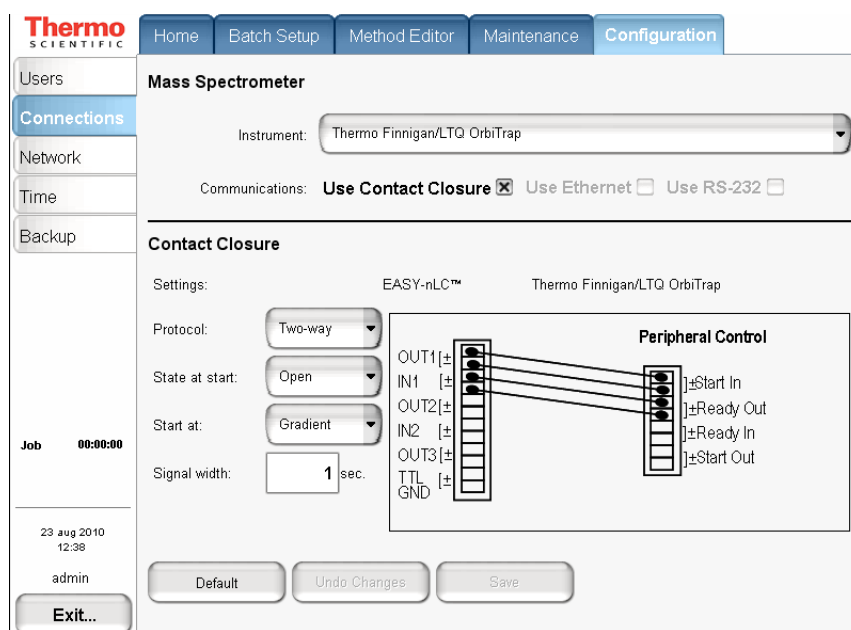
Setting Up the Mass Spectrometer Connection

IMPORTANT Before you connect the HPLC to a mass spectrometer, make sure that the protective grounding is shared between the instruments.

❖ To set up the mass spectrometer connection

1. Choose **Configuration > Connections**. The application displays a number of options.

Figure 21. Connections page



2. From the Instrument list, select the correct mass spectrometer (see [Figure 21](#)). The application comes preconfigured with most of the standard mass spectrometers used in proteomics, but if you cannot find your particular mass spectrometer, select Generic instead.

If your mass spectrometer does not appear on the list and you would like to see it supported in a future release of the EASY-nLC II application, provide this information to Thermo Fisher Scientific Customer Service (see [“Contacting Us”](#) on [page ix](#)).

3. Ensure that the **Use Contact Closure** option is selected for the HPLC-mass spectrometer communications protocol.
4. To define contact closure, from the Protocol list select **One-way** synchronization (see [Figure 21](#))—that is, a simple signal at the start of sample loading or start of gradient.

With a two-way connection, the HPLC waits for an acknowledgement that the mass spectrometer is ready to accept the sample. To correctly set the start state and signal width, refer to the mass spectrometer documentation for how to set up contact closure at the mass spectrometer end.

By setting the Start At value to be the sample loading start, you might be able to acquire data from early-eluting peptides that would not ordinarily bind to the column material and thus be missed from the actual gradient. For this to be effective, you must be running in a one-column system configuration mode.

Note Moving the Start At point to the sample loading start can lead to small retention time variations, due to possible differences in pressure buildup before the A/B mixing starts in the actual gradient.

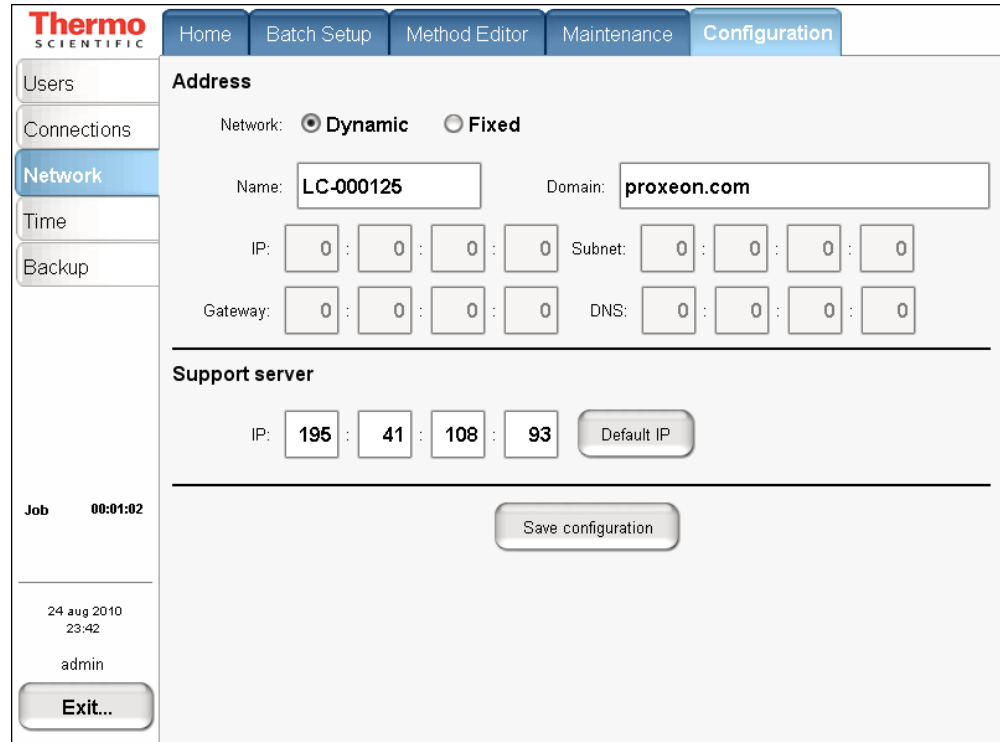
Network Connection

You can access the EASY-nLC II system from other computers (and for remote support) if it is connected to your network.

IMPORTANT Ask your local IT administrator to help you configure the EASY-nLC II system on the local network. The following information will help you through the process.

❖ **To connect the EASY-nLC II system to a local network**

1. From the top menu, press the **Configuration** tab and then press the **Network** tab. The Network view displays configuration address information.



2. For Network type, select **Dynamic** addressing (DHCP) or use a **Fixed** IP address. Your IT administrator knows which option to select.

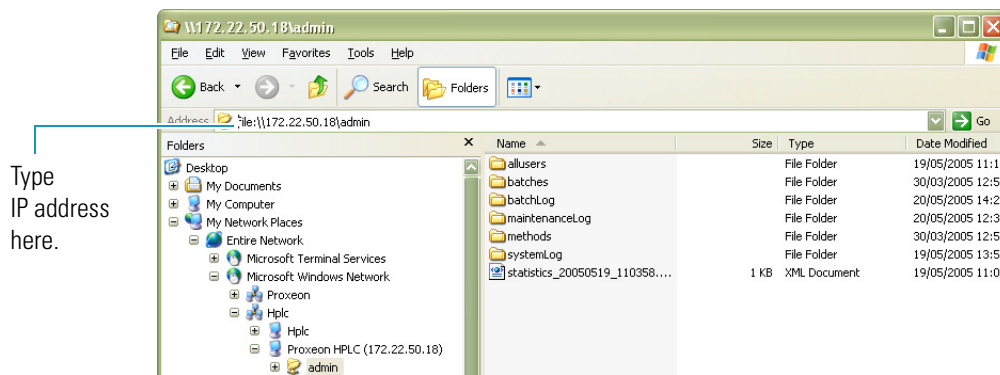
If IT asks about a MAC address for the instrument—that is, the instrument’s unique network identifier—it is labeled on the connector panel on the back of the instrument. The identifier is a series of numbers and letters and might look like this:

MAC address: 00:E0:4B:07:89:65

3. After setting up the network address, attach the Ethernet cable and restart the system to get connected. See [“Closing Down the EASY-nLC II System”](#) on page 33 for information about the power-down procedure, and then immediately go to [“Starting the EASY-nLC II System”](#) on page 19 and [“Logging In to the EASY-nLC II System”](#) on page 32.

External users can now access the file server in the EASY-nLC II system to look at log files, for example. Microsoft™ Windows™ Explorer might display the file server as shown in [Figure 22](#).

Figure 22. File server access through Windows Explorer



4. Type the IP address of the instrument, for example, **file:\\172.22.50.18**, and press ENTER.

You need a valid user name and password to enter the file space. Your user name and password are the same as those on the EASY-nLC II system.

User Permissions

There are four types of users: Guest (default), Normal, Super user, and Administrator.

Each user type has access to different levels of functionality in the EASY-nLC II application.

Table 8. User types and permission levels

Type/Menu	Guest	Normal	Super user	Administrator
Home	(X) ²	X ¹	X	X
Batch Setup	(X)	X	X	X
Method Editor	(X)	X	X	X
Maintenance			X	X
Configuration				X

¹ X = User has read-write access.

² (X) = User has read-only access.

Blank = Feature is not accessible.

Having a Guest user account means you can access the top menu but not enter data or issue commands to the system.

If a feature is not available, the application grays (blanks) it out, so the user cannot access that function.

Changing the Administrator Password

The EASY-nLC II application supplies a simple administrator password. Change it to a more secure password as soon as possible.

❖ To change the administrator password

1. Log in as **admin** (otherwise, see “Logging In to the EASY-nLC II System” on page 32).
2. In the top menu, press the **Configuration** tab and then press the **Users** tab. The Users page opens (Figure 23).

Figure 23. Users page under Configuration

The screenshot shows the Thermo Scientific EASY-nLC II web interface. At the top, there is a navigation bar with tabs: Home, Batch Setup, Method Editor, Maintenance, and Configuration. The Configuration tab is selected. On the left, there is a sidebar with tabs: Users, Connections, Network, Time, and Backup. The Users tab is selected. The main content area displays the configuration for the 'admin' user. The 'User' field contains 'admin'. The 'Name' field contains 'Administrator'. The 'Password' and 'Re enter' fields are masked with asterisks. The 'Level' is set to 'Admin' (radio button selected). Below the level, there are four dropdown menus for 'Graphs': Pump A, Pump B, Gradient, and Sample Cooling. At the bottom, there are buttons for 'Open', 'Close', 'Save', 'Delete', and 'New'. In the bottom left corner, there is a status bar showing 'Job 00:00:00', the date '24 aug 2010 17:17', the user 'admin', and an 'Exit...' button.

3. Press **Open** and select **Admin** from the User list. You can now enter and confirm a new password.
4. Press **Save** to save the user account information.

Creating New User Accounts

Use the Users page to create new user accounts for all the people or groups that will use the instrument. Each user has a private file space for batch jobs and methods that other users (apart from system administrators) cannot see.

In addition, there is a public file space for methods and batch jobs that all users can see and use. Only Super users and Administrators can copy files into the public spaces and only Administrators can delete files and directories from the public space.

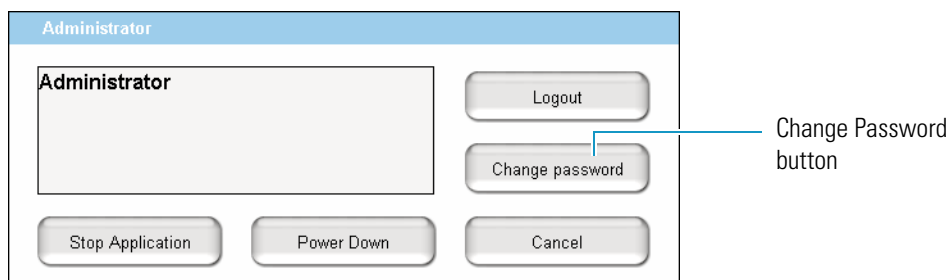
❖ To create a new user account

1. Log in as **admin** (otherwise, see “[Logging In to the EASY-nLC II System](#)” on [page 32](#)).
2. Choose **Configuration > Users**. The Users page opens. See [Figure 23](#) on [page 39](#).
3. Press **New**. Select which privileges to give the new user. For user privileges by user type, see [Table 8](#) on [page 38](#).

If you want the user to carry out instrument maintenance, you must select Super user as the user type. If you want the user to restore data to the system, select Administrator as the user type.

4. Create an initial password for the defined user.

Tip Only certain users have access to the Configuration > Users screen to change their own password. All users, however, can change their current password from the Logout confirmation box by pressing Change Password.



Configuring Column Setup, Loop Volume, Idle Flow, and IFC

Use the Devices page on the Maintenance tab to set global instrument configuration settings such as one- or two-column setup, loop size, idle flow settings, and IFC. For parameter descriptions, see [Table 9](#).

❖ To configure devices

1. Choose **Maintenance > Devices**.
2. From the Devices list, select **EASY-nLC II (HPLC)**.

3. Press the **Properties** tab to configure your settings (Figure 24).

Figure 24. Device configuration for HPLC on the Properties tab

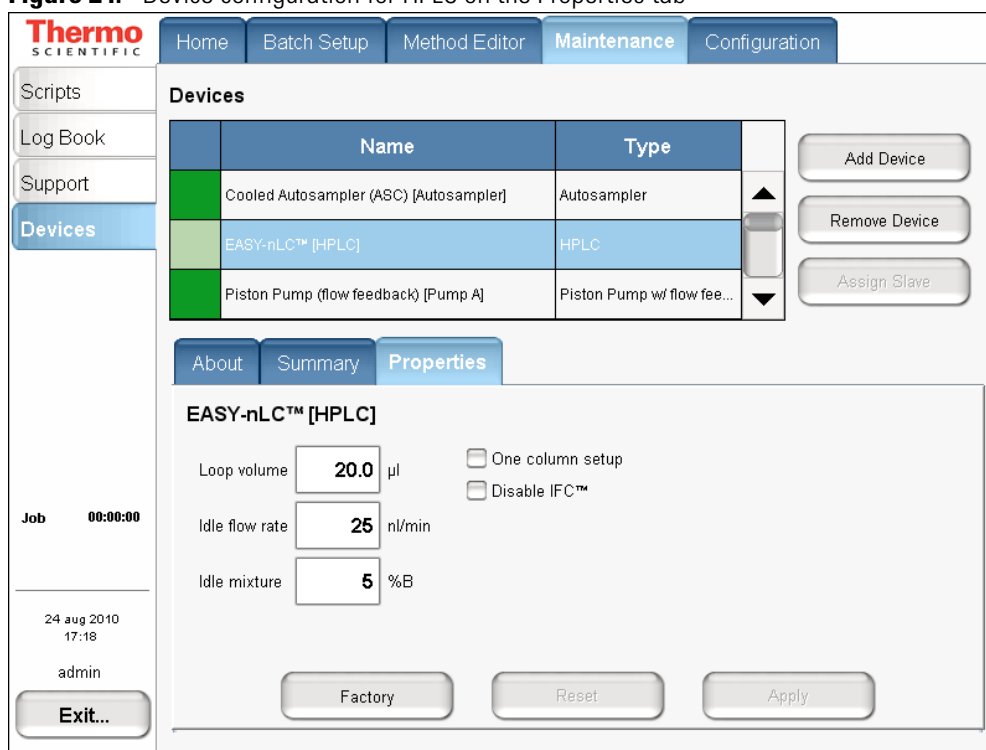


Table 9. HPLC parameter descriptions (Sheet 1 of 2)

Parameter	Description
Loop Volume	Make sure to configure the loop volume correctly, as method input parameters are limited by the loop volume. If a smaller loop is physically mounted over what is configured in the EASY-nLC II application, the instrument might actually draw sample into the pump.
Idle Flow Rate	Maintaining long-term emitter stability (with glass emitters) might require constant flow through the system. Idle flow is a functionality that activates when the last sample in a batch is done and is at a constant flow through the system at a specified rate and as a percentage of B. If you do not want such a flow, type zero (0) in the Idle Flow Rate box.

Table 9. HPLC parameter descriptions (Sheet 2 of 2)

Parameter	Description
One Column Setup	You can configure the EASY-nLC II instrument as a one-column or a two-column setup. If the system is operated as a one-column system, you can select its corresponding check box here, and the diagram on the Overview page of the Home menu will reflect this. The actual execution of the method is unaffected by this setting.
Disable IFC	For a description of IFC, see “Pump Flow Control” on page 5 . You can disable it by selecting this check box.

Preparation for Use

This chapter describes how to prepare the EASY-nLC II instrument for use.

Log in to the system with either Super user or Administrator privileges. For details, see “Logging In to the EASY-nLC II System” on page 32.

Contents

- [Preparing Solvent Bottles and Waste Containers](#)
- [Executing Maintenance Scripts](#)
- [Purging and Flushing the Pumps](#)
- [Installing Columns](#)

Preparing Solvent Bottles and Waste Containers

Before shipment, the EASY-nLC II system is flushed with methanol.



CAUTION Methanol (CAS number: 67-56-1) is highly flammable, and toxic by inhalation, ingestion, or skin absorption. Take appropriate measures to protect yourself and your equipment. Make sure that the mobile phases are miscible with methanol, or start up with an intermediate solvent as mobile phase.

Solvent A and B Bottles

Follow these recommended steps for preparing the two solvent bottles.

❖ To prepare the solvent bottles

1. Fill the 25 ml Schott Duran™ blue cap bottles with the mobile phase and degas them either by sonification or by sparging them with helium gas.
2. Place the bottles in the holder on top of the instrument: solvent A bottle in the front and solvent B bottle in the back position.
3. Mount the inline filter assembly on the blue cap lids as shown in [Figure 25](#) and [Figure 26](#):

Figure 25. Inline filter assembly (exploded view)

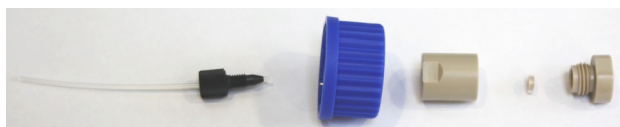
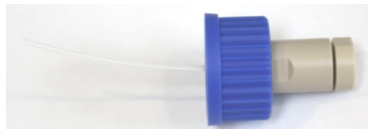


Figure 26. Assembled inline filter



IMPORTANT To establish proper surface wetting, prime the filters with methanol or acetonitrile.

Autosampler Bottles

The EASY-nLC II autosampler holds up to four bottles: three for washing solvents and the fourth for ejecting waste and cleaning the outside of the injection needle. For the bottle holder positions, see [Figure 27](#) on [page 45](#).

You can operate the instrument after installing the washing bottle for needle cleaning (position W4) and the solvent A bottle (position W3).

❖ To prepare the autosampler bottles

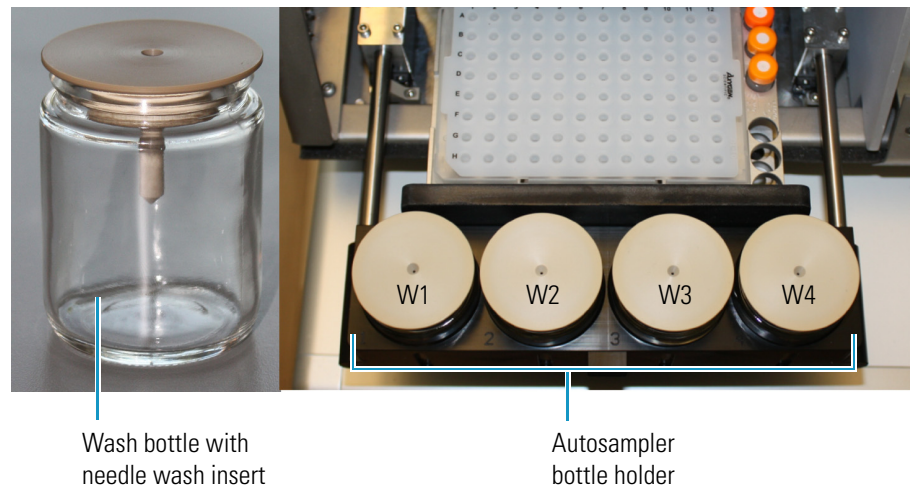
1. Fill one of the bottles with solvent A and degas it.
2. Turn on the system, go to **Home > Overview**, and press **Eject Plate**.
3. Put the bottle containing the needle wash insert into position W4.



CAUTION You must place the washing bottle holding the needle wash insert in position W4. Any other position could lead to instrument malfunction and possible breakdown.

- Put another bottle containing solvent A into position W3. The S pump is filled from this position. You can place wash solvents for customized washing procedures in position W1 and W2. See [Figure 27](#) on [page 45](#).
- Make sure all the bottles are installed with the lids fitting tightly.
- After positioning the bottles, press **Insert Plate** to move the plate back into the instrument.

Figure 27. Wash bottle with needle wash insert and view of the four positions on the autosampler bottle holder (ASC model)



Waste Container

The EASY-nLC II waste container sits behind a small door next to the autosampler door. All solvent waste, except waste from the autosampler, goes into this waste container.

❖ To prepare the waste container

- Open the autosampler door and insert the waste container in the compartment (see [Figure 28](#)), without mounting the yellow lid.

Figure 28. Waste container compartment



- Make sure to properly position the waste container.
- Close the door.

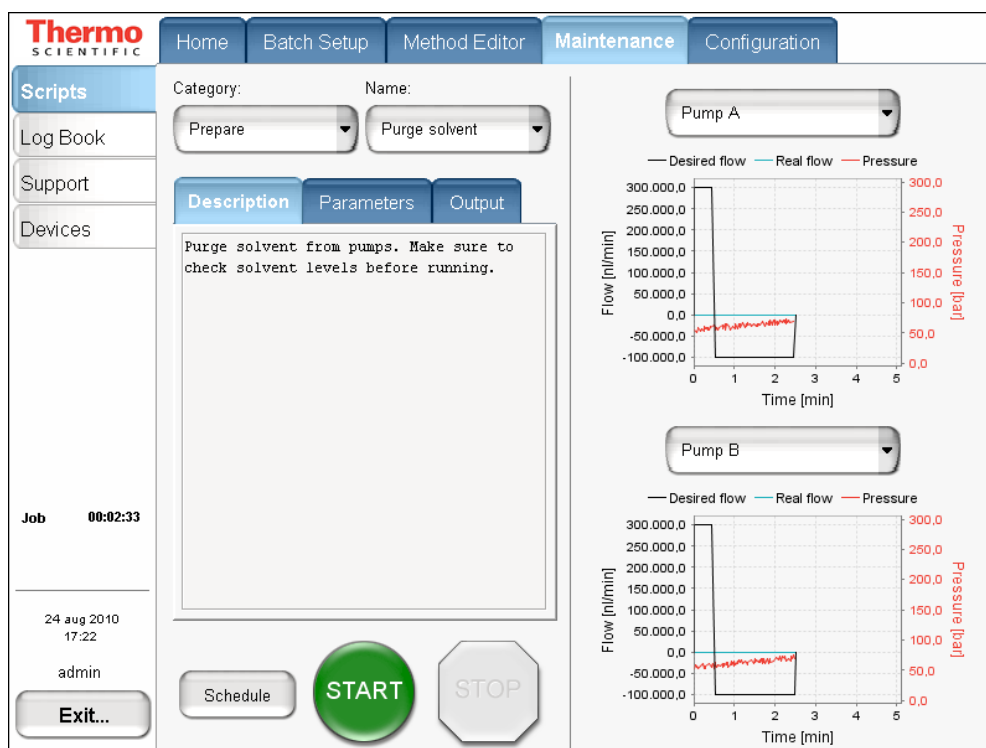
Executing Maintenance Scripts

Note

Before proceeding with the preparation for use, make sure to perform all tasks described so far in “Preparing Solvent Bottles and Waste Containers” on page 43.

You cannot execute maintenance scripts when the EASY-nLC II instrument is running batches.

Figure 29. Scripts page under Maintenance



The green START button is active when a job or script is ready to execute. During execution you can terminate a running job by pressing the now active (red) STOP button.

You can automate the start of certain scripts, for example, scheduling them for later execution. However, you can schedule only those scripts where the Schedule button is active. You cannot schedule some scripts because they require supervision or hardware changes to the system.

Purging and Flushing the Pumps

Because the pumps have been purged with methanol before delivery of the EASY-nLC II instrument, make sure to purge them thoroughly before doing any analytical runs. In addition, you must purge the pumps when switching to a new solvent.

The purge pump job fills the chosen pump or pumps with solvent and afterwards ejects it into the waste container. You can specify how many iterations of this job you want performed.

After purging the pumps, flush them to remove any air that might be trapped inside. You can specify the maximum number of iterations for this job, the period of time that the pumps are being pressurized, and the success criteria for stopping the operation earlier than scheduled.

❖ To purge and flush pumps A and B

1. Go to **Maintenance > Scripts** to open the Scripts view.
2. In the Category list, select **Prepare**, and in the Name list, select **Purge Solvent**.

The screenshot displays the 'Scripts' view in the Thermo Scientific software. The 'Category' is set to 'Prepare' and the 'Name' is 'Purge solvent'. The 'Parameters' tab is selected, showing a table with the following data:

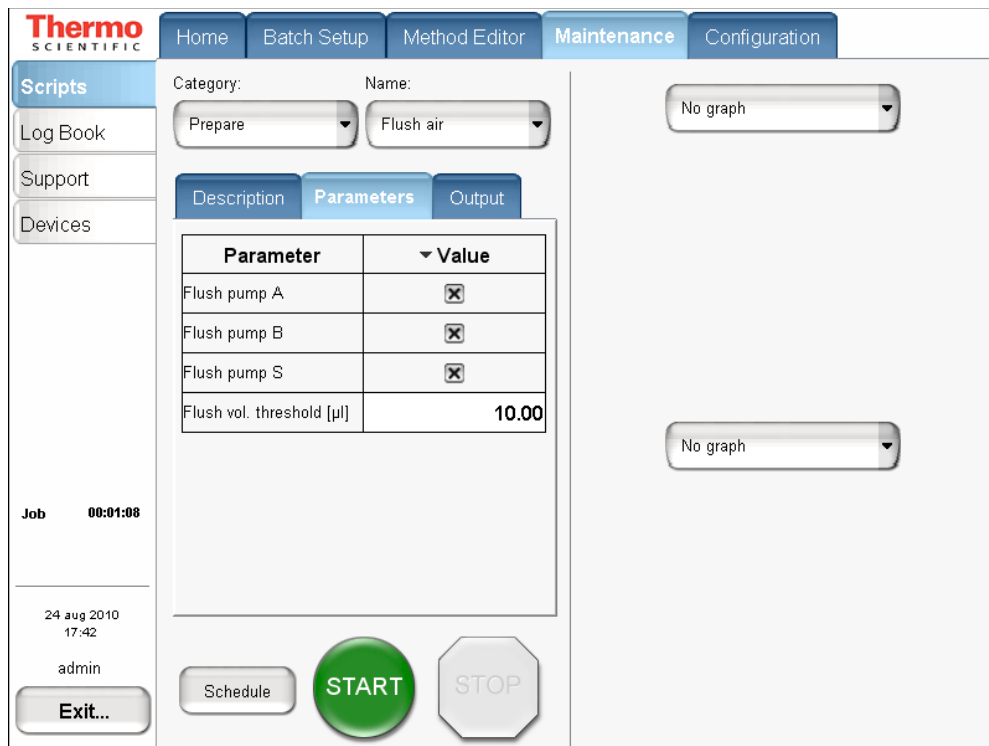
Parameter	Value
Purge iterations	10
Purge pump A	<input checked="" type="checkbox"/>
Purge pump B	<input checked="" type="checkbox"/>
Purge pump S	<input checked="" type="checkbox"/>

Below the table are buttons for 'Schedule', 'START', and 'STOP'. On the right, there are two graphs for 'Pump A' and 'Pump B'. Each graph plots 'Flow (ml/min)' on the left y-axis (0.0 to 300.000.0) and 'Pressure (bar)' on the right y-axis (0.0 to 300.0) against 'Time (min)' on the x-axis (0 to 5). The graphs show 'Desired flow' (black line), 'Real flow' (blue line), and 'Pressure' (red line).

3. Press the **Parameters** tab, select the cell next to Purge Iterations, and type **10**. This is the number of times that the pumps will completely refill and eject their volume.

Note Bottle three (W3) in the autosampler bottle holder fills the S pump, and the solvent ejects into bottle four (W4), which is always used for waste liquid.

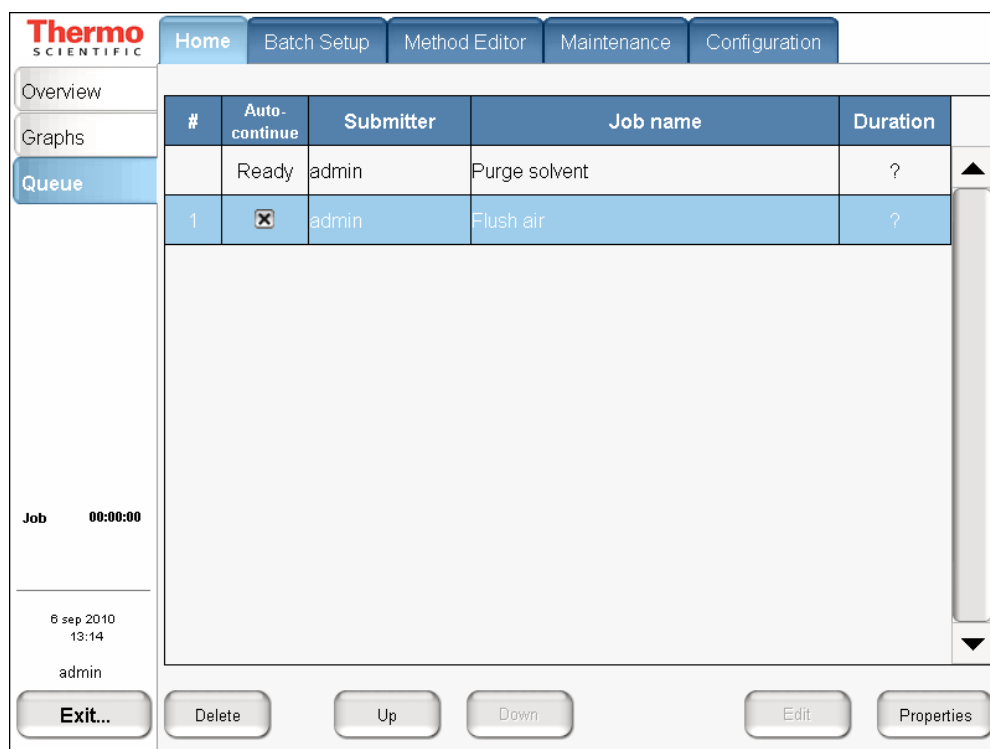
4. To save this setting and continue, press **Schedule**. The Home > Queue page appears.
5. Return to **Maintenance > Scripts**.
6. In the Category list, select **Prepare**, and in the Name list, select **Flush Air**.



7. Press the **Parameters** tab, select the cell next to Flush Vol. Threshold [µl], and type **10**. This is the acceptable amount of compressibility that helps ensure good system performance, overall.

8. Press **Schedule**.

The Home > Queue page appears.



9. In the Jobs list, make sure the check box in the Auto-Continue column is selected for **Flush Air**.

10. Choose **Home > Overview** again and press **START**.

Default graphs are selected in the two graph windows in Maintenance > Scripts, but you may change the selection from the corresponding lists. These graphs help you to monitor the job.

11. Wait for the job to end (the flush air script continues until the flush air volume values are below the specified threshold).

Note While a maintenance job or a script is running, you cannot start other jobs or scripts.

Installing Columns

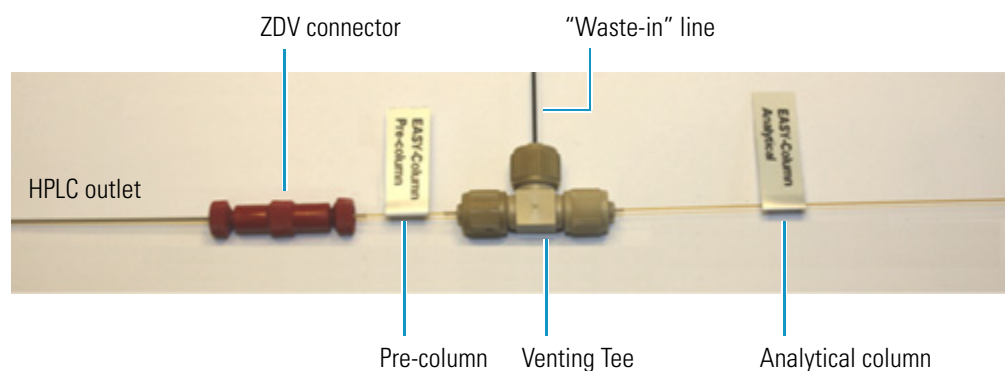
The EASY-nLC II instrument comes with a trap column (EASY-Column™, 2 cm, ID 100 µm, 5µm, C18-A1), an analytical column (EASY-column, 10 cm, ID 75 µm, 3 µm, C18-A2), and the unions and fittings for connecting 375 µm OD and 1/32 in. OD columns. You can use other types of column ODs. Contact Thermo Fisher Scientific for help choosing the appropriate fittings.

Two 1/32 in. OD PEEKsil™ lines are attached to the instrument, one called “col out” and the other, “waste in.”

❖ To connect columns

1. Connect the red union (SC600) on the “col out” PEEKsil. Read the “Instructions for Use” note provided with the union.
2. Connect the pre-column to the other end of the union, using the provided sleeves (SC603) if it is a 375 µm OD column.
3. Connect the flow-out end of the pre-column to the Venting Tee (SC601). Again, use the provided sleeve (SC603) if it is a 375 µm OD column.
4. Connect the angled end of the Venting Tee with the “waste in” PEEKsil.
5. Connect the straight end of the Venting Tee to the analytical column and use the provided sleeve (SC603) if it is a 375 µm OD column.
6. Connect the analytical column outlet to the transfer line going to the detector.

Figure 30. EASY-nLC II column setup (375 µm OD columns)



Note The systems maximum pressure is set at the factory to 300 bar (280 bar with IFC enabled). Parameters such as column resin and frit size can vary greatly and generate different back pressure.

You can perform reequilibration and sample loading quicker at higher flow rates, which help to reduce cycle time when analyzing samples.

Preparing the Pre-Column

This procedure assumes that IFC is activated (the default). If IFC is disabled, the EASY-nLC II application ignores the Max Pressure input field and, instead, the set flow passively regulates the pump.

❖ To prepare the pre-column

1. Choose **Maintenance > Scripts**.
2. Select **Prepare** in the Category list.
3. Select **Precolumn equilibration** in the Name list.

The screenshot displays the Thermo Scientific software interface. At the top, there are navigation tabs: Home, Batch Setup, Method Editor, Maintenance (selected), and Configuration. On the left, a sidebar contains 'Scripts', Log Book, Support, and Devices. The main area is divided into 'Category' (Prepare) and 'Name' (Precolumn equilibration). Below this, there are three tabs: Description, Parameters (selected), and Output. The Parameters tab contains a table with the following data:

Parameter	Value
Volume [µl]	30.00
Flow [µl/min]	5.00
Max pressure [bar]	250.00

Below the table are buttons for 'Schedule', 'START' (highlighted in green), and 'STOP'. On the right side, there are two graph windows for 'Pump A' and 'Pump B'. Each graph plots 'Flow [ml/min]' (left y-axis, 0.0 to 300,000.0) and 'Pressure [bar]' (right y-axis, 0.0 to 300.0) against 'Time [min]' (x-axis, 0 to 5). The graphs show 'Desired flow' (black line), 'Real flow' (blue line), and 'Pressure' (red line).

4. In the Scripts view, press the **Parameters** tab and set the desired flow, volume, and maximum pressure. Thermo Fisher Scientific recommends running the reequilibration at a set pressure rather than a set flow, to make sure the column pressure limit and the instrument pressure limit (300 bar) are not exceeded. If the flow field is left empty, the pump will run at the specified pressure. For best results, use a reequilibration volume of 10 column volumes.
5. Press **START**. The default graphs appear in the two graph windows, but you may change the selections from their corresponding lists. These graphs help you to monitor the progress.
6. Monitor the actual flow rate to become familiar with the flow/pressure relationship on your particular column.

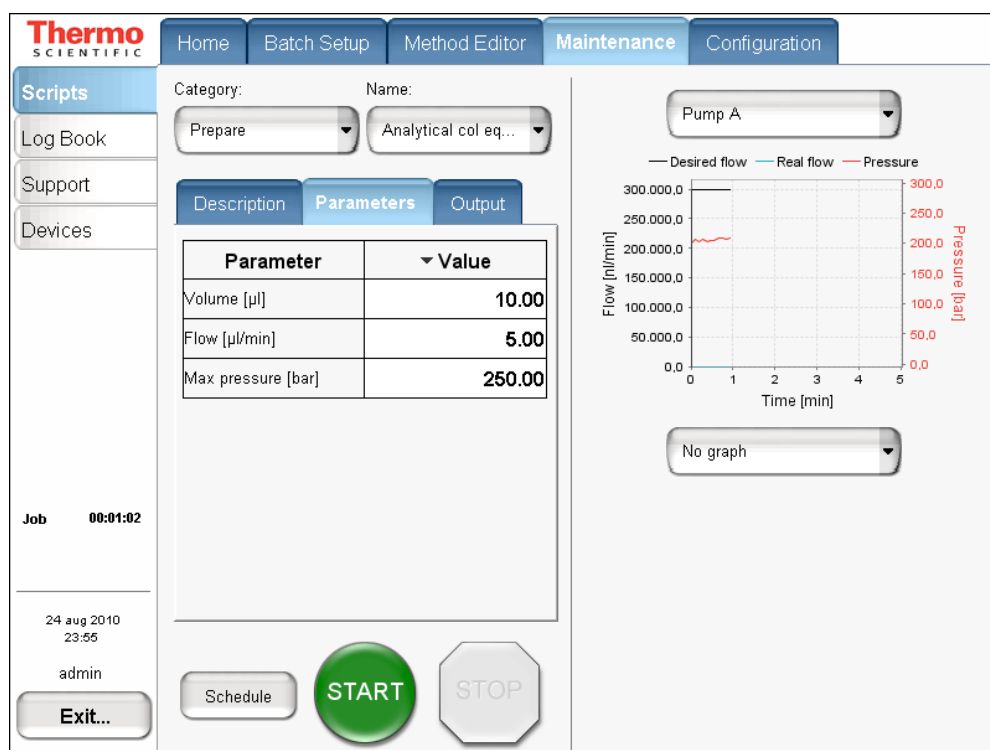
Equilibrating the Analytical Column

This procedure assumes that IFC is activated (the default). If IFC is disabled, the EASY-nLC II application ignores the Max Pressure input field and, instead, the pump is regulated passively by the set flow.

❖ To equilibrate the analytical column

1. Choose **Maintenance > Scripts** (see Figure 29).
2. Select **Prepare** in the Category list.
3. Select **Analytical col equilibration** in the Name list.

Figure 31. Scripts page under Maintenance



4. In the Scripts view, press the **Parameters** tab and set the desired flow, volume, and maximum pressure. Thermo Fisher Scientific recommends running the equilibration at a set pressure rather than a set flow, to make sure the column pressure limit and the instrument pressure limit (300 bar) are not exceeded. If the flow field is left empty, the pump runs at the specified pressure. For best results, use an equilibration volume of 10 column volumes.
5. Press **START**. The default graphs appear in the two graph windows, but you may change the selections from their corresponding lists. These graphs help you to monitor the progress.
6. Monitor the actual flow rate to become familiar with the flow/pressure relationship on your particular column or columns.

Running the First Sample

This chapter describes how to perform analysis using the EASY-nLC II system. Thermo Fisher Scientific recommends that you quickly start using the application to run additional analytical tests, as practical experience is the most efficient way to understand an instrument such as an HPLC.

Contents

- [Preparing a Well Plate](#)
- [Creating a Method](#)
- [Creating a Batch](#)
- [Starting Sample Acquisition](#)
- [Monitoring the Run](#)
- [Stopping Sample Acquisition](#)
- [Editing the Running Batch](#)
- [Troubleshooting a Sample Run](#)

Preparing a Well Plate

If you have not made a plate selection, go to “[To select a given plate format](#)” on page 76.

The EASY-nLC II instrument ships with adaptors for vials (56 vials in a 6×8+6 configuration), MTP (96- and 384-well), and PCR strips.

Note The instrument comes preinstalled with a 20 µL loop. With this loop size, the instrument is capable of picking up sample volumes between 0 and 18 µL.

❖ To prepare the well plate

1. Load some vials/wells with your sample, preferably a known standard.
2. Make sure that no air is trapped in the sample vials/wells and that the sample is in the bottom of the vials/wells. To do this, tap the vial/plate gently against a hard surface to have the sample move to the bottom.
3. Put on the vial cap or plate mat.

Creating a Method

The Edit view under Method Editor consists of five pages to guide you through all the parameters required for a working LC method. The protocol is based on IFC being active (the default system). If you deactivate IFC, the EASY-nLC II application ignores the maximum pressure input fields and instead, the pumps are regulated passively by the set flow (only relevant for sample loading and equilibration steps).

The following sections describe the five pages and guidelines to creating a method.

1. Choose **Method Editor > Edit**.
2. If you are not already on page 1/5, switch pages by pressing the appropriate arrow near the bottom of the screen.

Page 1/5

❖ To open existing methods and close or save the currently opened method

1. Press **SAVE**.
2. In the Save File dialog box, press the white box beneath the name.
3. Type **method 1** on the keyboard and press **OK**.
4. Press **SAVE** in the Save File dialog box.

You may enter a detailed explanation or description in the white text box.

5. Go to page 2/5 by pressing the right arrow.

Page 2/5

Use page 2/5 (see Figure 32) to set the parameters for picking up the sample from the well plate and loading it onto the pre-column.

Figure 32. Page 2/5, Edit view in the Method Editor

The screenshot shows the Thermo Scientific Method Editor interface. At the top, there are navigation tabs: Home, Batch Setup, Method Editor (selected), Maintenance, and Configuration. Below the tabs, there is a 'File' menu and a 'Name' field containing 'BSA 30 min' with a 'Save' button to its right. The main area is divided into two sections: 'Sample pickup' and 'Sample loading'. In the 'Sample pickup' section, there are input fields for 'Volume: 10.00 µl' and 'Flow: 20.00 µl/min', accompanied by a schematic diagram of the autosampler. In the 'Sample loading' section, there are input fields for 'Volume: 30.00 µl', 'Flow: 7 µl/min', and 'Max. pressure: 250 Bar', accompanied by a schematic diagram of the sample loading process. On the left side, there is a 'Job' status indicator showing '00:00:07', a date and time '24 aug 2010 17:27', and the user 'admin'. An 'Exit...' button is located at the bottom left. At the bottom center, there are navigation arrows and the page number '2/5'.

❖ To set up the autosampler to pick up sample

1. In the Sample Pickup area, press the Volume box and enter the default amount of sample to be picked up. To prevent contamination of sample pump S, the maximum volume is set to loop size minus 2 µL.
2. Press the Flow box to set the flow rate at which the autosampler picks up sample.

The autosampler is able to pick up sample with a flow between 0 and 40 µL/min. Set the speed according to the sample viscosity and accuracy needed; however, 20 µL/min is a good starting point when sample is in an aqueous solution. No excess of sample will be picked up—just the exact amount of the requested sample.

❖ To set up sample loading

1. In the Sample Loading area, press the Volume box to set how much solvent A to use to load the sample from the loop onto the pre-column. Normally it is sufficient to load the pre-column with 2.5 times the amount of sample picked up.
2. Press the Flow box to set the flow rate for sample loading onto the pre-column. If you leave the input box empty, the pump operates at the set pressure (IFC enabled).

3. Press the Max. Pressure box to set the maximum allowed pressure. If you leave the input box empty, the pump operates at the set flow (as long as it is below the instrument pressure of 280 bar).

Note If you specify both a flow rate and a maximum pressure, the pump flow is limited by the parameter that is reached first.

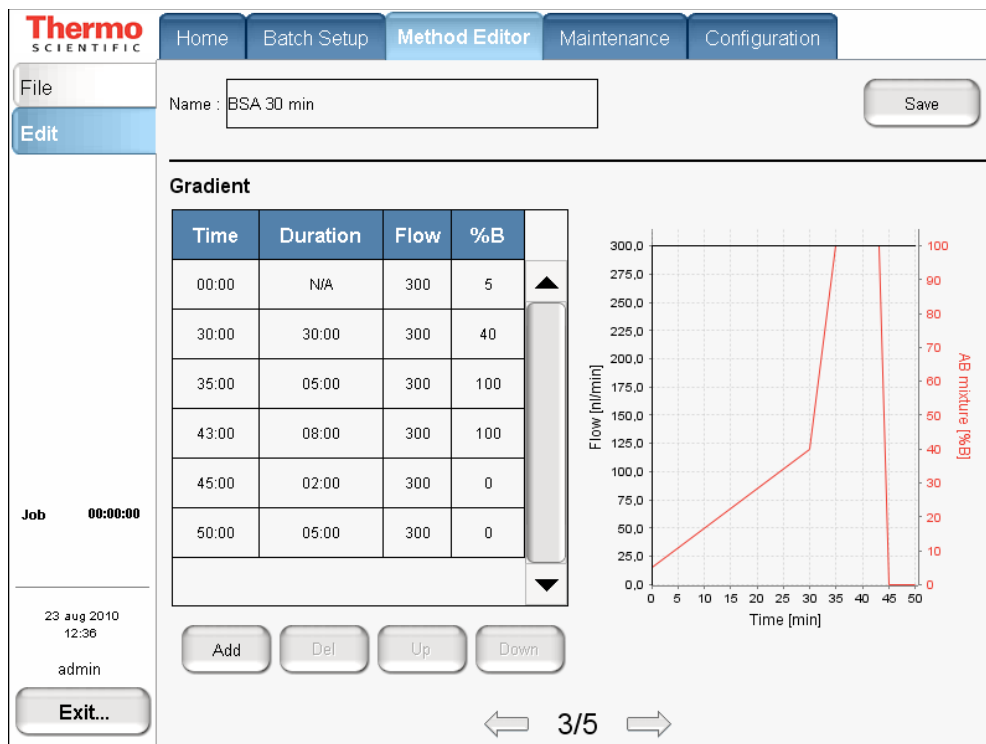
Page 3/5

Page 3/5 is for building a gradient. It consists of a schematic view for setting parameters and a graph view showing the gradient as a function of time (see [Figure 33](#)).

You can use the four buttons at the bottom of the page for adding lines to the gradient, moving lines up and down, and deleting lines. To move or delete lines, select the line by pressing it.

The rest of this section guides you through building a gradient for analyzing a BSA digest on a C18 column. Because this procedure is intended to help you set up a method, make sure to alter the parameters stated to match your own sample and column characteristics.

Figure 33. Page 3/5, Edit view in the Method Editor



❖ **To build a gradient**

1. Focusing on the first row of the gradient table, press the first cell in the Flow column, enter **300** nL/min in the pop-up box, and press **ACCEPT**.
2. Press the first cell in the % B column, enter **10** (%) in the pop-up box, and press **ACCEPT**.

The gradient is now set to start at 300 nL/min and 10 %B.

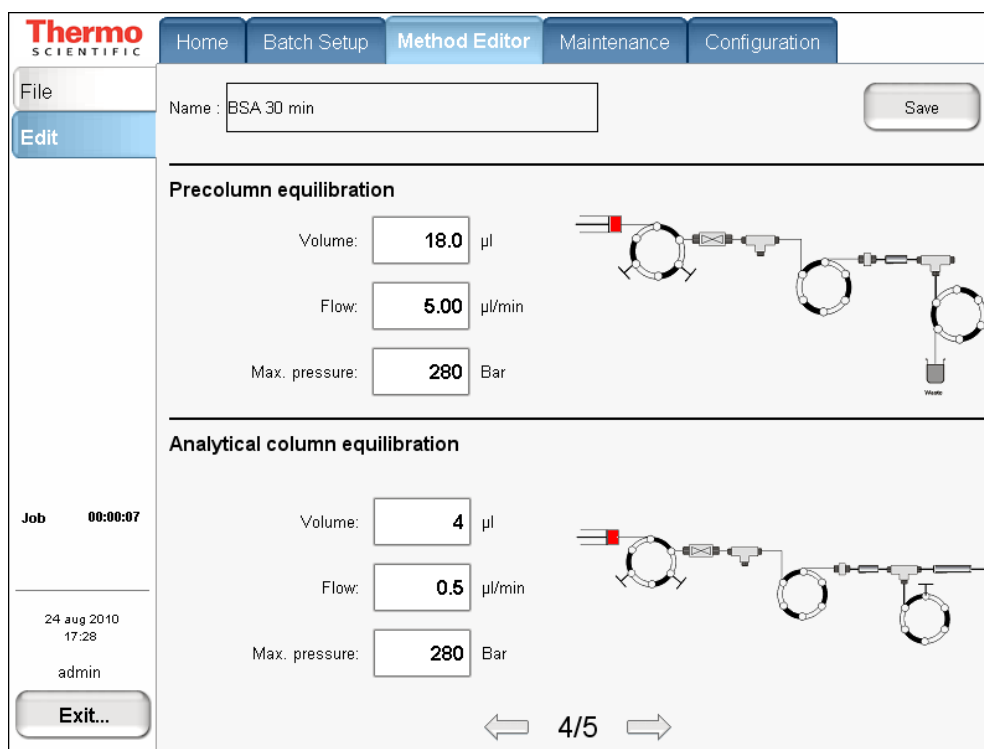
3. Press **ADD** to add another row.
4. Press the cell in the second row under the Duration column and set the duration to **20** min.
5. Press the cell in the second row under the Flow column and set it to **300** nL/min.
6. Press the cell in the second row under the % B column and set it to **40**.
7. Press **ADD** to add another row.
8. In this new row, set duration to **5** min, flow to **300** nL/min, and % B to **100**.
9. Press **ADD** to add another row.
10. In this new row, set the duration to **15** min, flow to **300** nL/min, and % B to **100**.

The gradient is now completed.

Note You do not need to reequilibrate the columns at the end of the gradient. The parameters for reequilibration are entered directly and separately as described in the next section. The reequilibration runs parallel with the sample pickup at the beginning of a sample execution run.

Use page 4/5 to set parameters for the column reequilibrations (see Figure 34).

Figure 34. Page 4/5, Edit view of the Method Editor



❖ **To set up the pre-column equilibration**

1. In the Precolumn Equilibration area, press the Volume box to set the amount of solvent A to use for reequilibration of the pre-column. For best results, use 10 column volumes for reequilibration.
2. Press the Flow box to set the flow rate for performing the reequilibration of the pre-column. If you leave the input box empty, the pump will operate at the set pressure.
3. Press the Max. Pressure box to set the maximum allowed pressure. If you leave the input box empty, the pump operates at the set flow (as long as it is below the instrument pressure of 280 bar).

Note If you specify both a flow rate and a maximum pressure, the pump flow is limited by the parameter that is reached first.

❖ **To equilibrate the analytical column**

1. In the Analytical Column Equilibration area, press the Volume box to set the amount of solvent A to use for equilibration of the analytical column. For best results, use 10 column volumes for reequilibration.

2. Press the Flow box to set the flow rate for performing the equilibration of the analytical column. If you leave the input box empty, the pump will operate at the set pressure.
3. Press the Max. Pressure box to set the maximum allowed pressure. If you leave the input box empty, the pump operates at the set flow (as long as it is below the instrument pressure of 280 bar).

Note If you specify both a flow rate and a maximum pressure, the pump flow is limited by the parameter that is reached first.

Page 5/5

Use Page 5/5 for setting up the autosampler washing procedures (see Figure 35). The EASY-nLC II instrument conducts the wash while the gradient is running.

Autosampler Wash

For most samples in low concentrations, flushing the needle and loop with solvent is sufficient. You can set the amount of solvent for this action in the Flush Volume box.

Use 100 µL as the default flush volume.

Note The minimum flush volume that you can set is the *loop volume + needle volume + 1 µL*.

Figure 35. Page 5/5, Edit view of the Method Editor

The screenshot shows the 'Method Editor' interface for 'Page 5/5'. The main area is titled 'Autosampler wash'. It features a 'Flush volume' input field containing '100.0 µL'. To the right is a schematic diagram of the autosampler. Below the flush volume is a 'Custom wash' section with an unchecked checkbox. Underneath is a table with the following structure:

Order	Source	Volume	Cycles

At the bottom of the interface, there are five buttons: 'Add line', 'Delete line', 'Copy line', 'Up', and 'Down'. The page number '5/5' is centered at the bottom.

Custom Wash

For a more thorough wash of the needle and loop, you can select the Custom Wash check box. Use this option to set up a user-definable procedure for washing the loop and needle.

When you use a custom wash, the flush volume that you set in the Autosampler Wash area defaults to a preset volume to flush the loop and needle with as much solvent as possible before starting the custom wash.

The principle for setting the Custom Wash parameters is similar to that for setting up the gradient. The Custom Wash table consists of three user-editable parameters described in [Table 10](#).

Table 10. Custom wash parameters

Parameter	Description
Source	Defines which wash bottle is picking up the solvent. You can choose between position 1, 2, or 3 (position 4 is for waste and cleaning of the outside of the needle). Position 3 is the default position where the pump is filled and normally contains solvent A. For example, you can fill position 1 and 2 with other solvents and select them as the Source to better clean the needle and loop.
Volume	Defines how much solvent is picked up. To avoid any contamination between the solvent being picked up and the pump, the pump is set to pick up 1 μ L of air before picking up sample, and the volume maximum is set to the following: $\text{needle volume} + \text{loop volume} - 2 \mu\text{L}$
Cycles	Defines how many cycles are going to be performed. Every cycle is completed with the volume set under volume being flushed out in wash bottle 4.

When the custom wash has ended, the autosampler automatically empties the pump and refills with solvent from bottle 3 to be ready for the next injection.

Creating a Batch

Use the Edit view under Batch Setup to set up and schedule batches for sample acquisition on the EASY-nLC II system. (See also “[Preparing a Well Plate](#)” on [page 54](#) and “[Creating a Method](#)” on [page 54](#).)

❖ To create a batch with the Microtiter plate

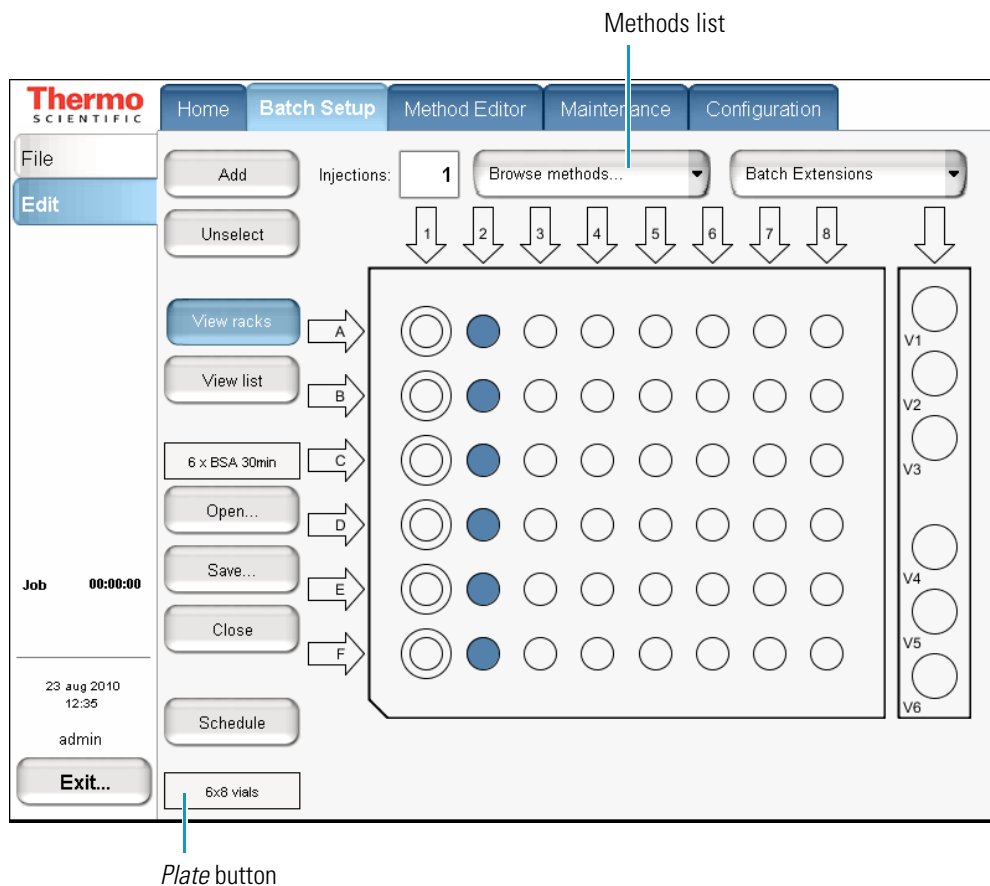
1. Go to **Batch Setup > Edit**. If you do not see a plate overview, press **View Racks** (see [Figure 36](#) on [page 61](#)).

2. Browse the methods list next to the Injections box, and select **method1** (created in “Creating a Method” on page 54).
3. If the plate format is not set to your installed plate, choose an appropriate plate by pressing the *plate* button. The *plate* button is the rectangular button imprinted with the plate name in the lower left side. The default plate is 6 × μl8 vials.
4. Press the DOWN ARROW labeled “1” to select all vials/wells from the first column (A1 to H1).

Note You can also select vials/wells by pressing them individually. You can clear well selections by pressing them again.

5. Press **Add** to link the selected vials/wells with the chosen method. Your touch-screen should look similar to the screen in Figure 36.

Figure 36. Vial/Well selection in the Edit view under Batch Setup



Note To use different methods in the same batch, choose another method, select new wells, and press **Add**.

6. Press **Save** to save the batch.

7. Press **Schedule** to submit the batch to the queue.

The application switches to the Home > Queue page (see [Figure 37](#)).

Figure 37. Queue page under the Home tab

#	Auto-continue	Submitter	Job name	Duration
	Ready	admin	6 x BSA 30min	11:04:15
1	<input type="checkbox"/>	admin	admin-20100824-1714	11:04:15

8. If necessary, move any scheduled batches in the job queue by selecting it and pressing **Up** or **Down**. You can also delete batches from the Queue by selecting the batches and pressing Delete.

For a batch run to start automatically after the previous batch is finished, you must select its corresponding check box in the Auto-Continue column. In the case in [Figure 37](#), the admin-20100824-1714 job does not auto-continue when the 6 x BSA 30min job is done.

Starting Sample Acquisition

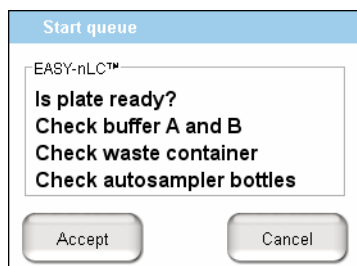
Before starting the sample acquisition, make sure that the system is properly set up. This includes having contact closure (see [“Connecting to the Mass Spectrometer through Contact Closure”](#) on page 13). Also make sure that the mass spectrometer or any other detector has been set up in accordance with the method.

At this point, you might want to run the maintenance script outlined in [“Test – MS Connection”](#) on page 111.

❖ To verify that the mass spectrometer connection (contact closure) works as expected

1. Choose **Home > Overview**.
2. Press **START** to start running the batch. A message box opens, prompting you to check that various conditions are fulfilled (see [Figure 38](#)).

Figure 38. Start Queue message box



3. Press **Accept**. The instrument now starts processing the batch.

Monitoring the Run

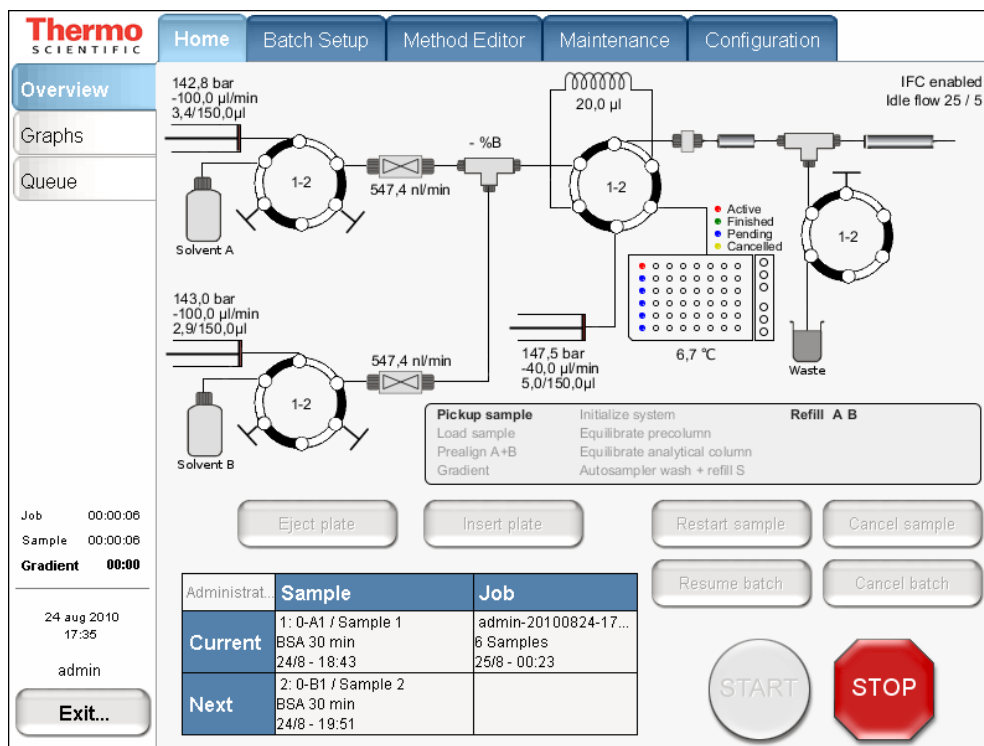
You can monitor the run and the system status during analysis from two different screens: Home > Overview or Home > Graphs.

- [Using the Overview Page in the Home Menu](#)
- [Using the Graphs Page in the Home Menu](#)

Using the Overview Page in the Home Menu

The Overview page provides a full schematic overview of the system, complete with real-time updating of all critical component data (Figure 39).

Figure 39. Overview page under Home



- The system schematic provides information on all hardware parts in the system:
 - Pump: position, pressure, and flow
 - Valve: position
 - Percentage of B being delivered
 - Actual flow being measured
 - Samples to be analyzed, plus their current status
 - Current sample and job, plus the next sample and job
- The sample analysis process can be divided into different activities such as Pickup sample, Load precolumn, and so on. When executed, these activities are highlighted in the activity box in the center of the page and give you a quick overview of the sample's progress.
- When the batch starts, the green START button is grayed out and the red STOP button becomes active.
- When the instrument is running, three counters measure the time for job, sample, and gradient.

Using the Graphs Page in the Home Menu

The Graphs page displays up to four graphs with historic data for the currently active sample (Figure 40).

Figure 40. Graphs Page under Home

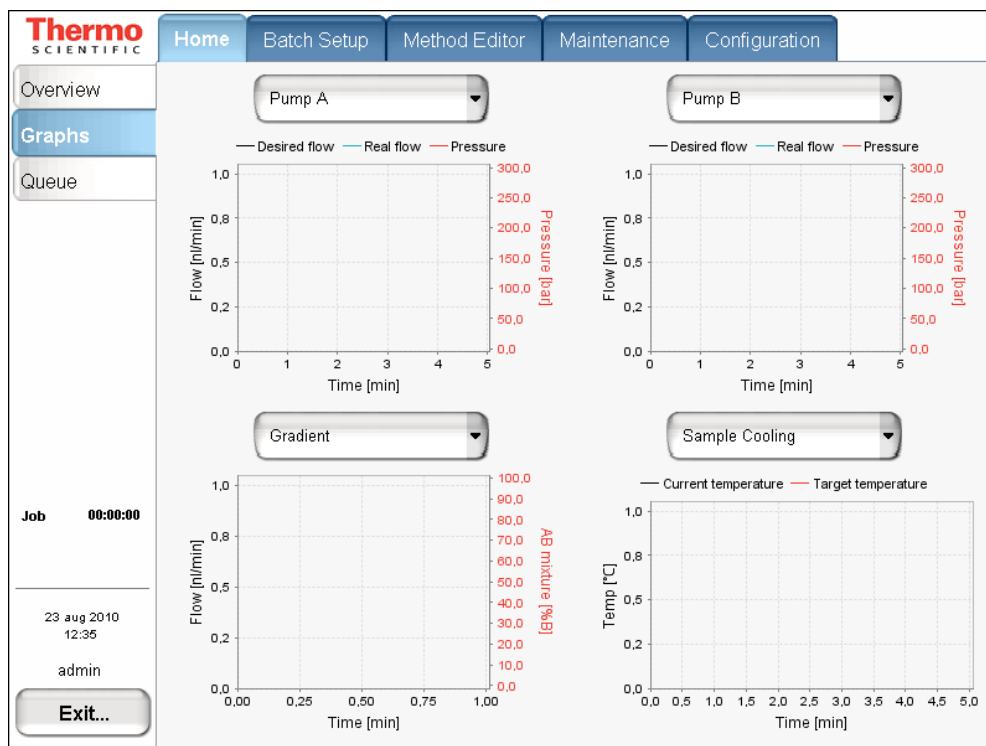


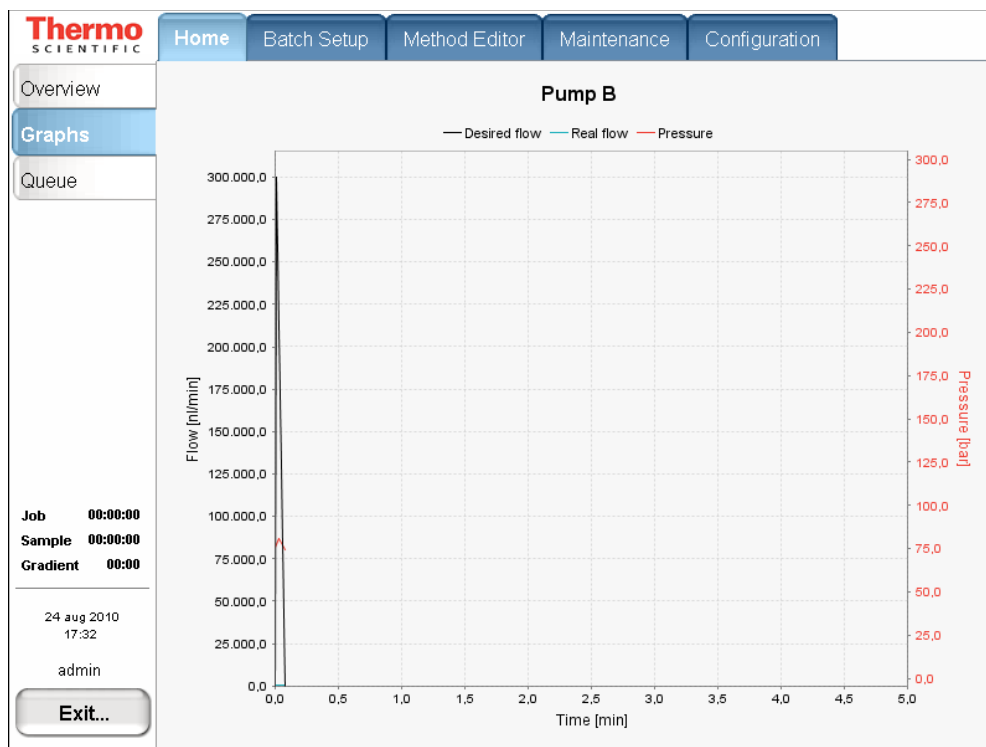
Table 11 describes the different types of graphs that are available. To change graphs, select from their corresponding lists.

Table 11. Graphical data types

Graph	Description
Pump (3)	For each of the three pumps, shows desired flow (black line), real flow (blue line), and pressure (red line) over time.
Valve (4)	For each of the four valves, shows valve position over time.
Gradient	Shows the theoretical gradient (dotted red line) and the actual gradient dynamically calculated on the feedback values from the pumps (full red line), plus the actual flow calculated on the feedback values from the pumps.
Temperature	Shows the measured temperature on the plate cooler over time.

To see a more detailed view, you can touch the graphs to enlarge them, for example, the flow graph for Pump B in Figure 41. Touch them again to minimize the view. Also, by touching the *y* axis you can cycle through several preset scales.

Figure 41. Enlarged graph on the Graphs page



Stopping Sample Acquisition

You can stop the sample acquisition during the run by pressing the red STOP button on the Home > Overview page. This halts all actions on the HPLC system and the following buttons become active:

- **Restart Sample:** Runs the method from its beginning again, cleaning up the system as it does. Only press this button if you are sure you have enough sample material in the Microtiter plate.
- **Cancel Sample:** Cancels the current sample and skips to the next sample in the batch job.
- **Cancel Batch:** Cancels the entire current batch job.
- **Eject Plate:** Ejects the plate for your inspection to help you determine how to proceed.

Editing the Running Batch

You can edit the running batch if you want to add sample or remove samples that are not yet tested.

❖ **To edit the running batch**

1. Choose **Home > Queue**.

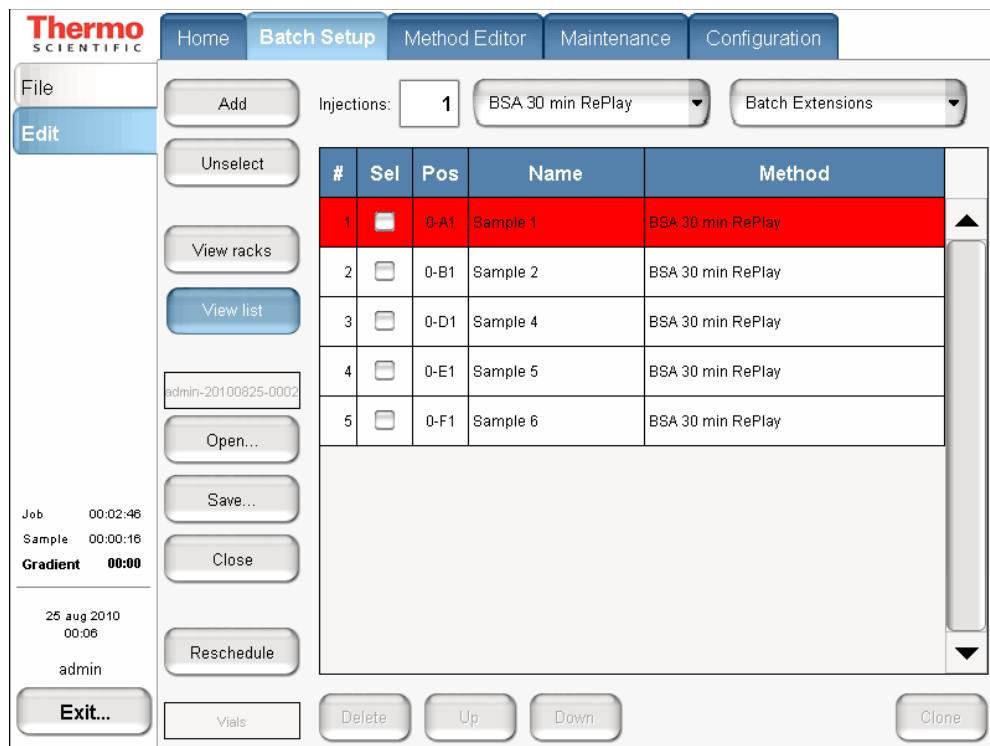
#	Auto-continue	Submitter	Job name	Duration
1	Running	admin	admin-20100825-0002	06:55:00
1	<input type="checkbox"/>	admin	6 x BSA 30min	06:48:00

Job 00:02:38
 Sample 00:00:08
Gradient 00:00
 25 aug 2010 00:05
 admin

Exit... Delete Up Down Edit Properties

2. Select the batch job from the list and press **Edit**. The Edit page under the Batch Setup tab appears (see Figure 42).

Figure 42. Edit page under Batch Setup tab



3. Edit the batch—that is, add, change, or remove samples that are not yet tested.
4. Press **Reschedule**. The batch continues running.

5. (Optional) Go to **Home > Queue** (see Figure 43), check the sample list by pressing **Properties**, and then press **Show Samples**.

Figure 43. Sample list on the Queue page of the Home tab

The screenshot displays the Thermo Scientific software interface. At the top, there is a navigation bar with tabs: Home, Batch Setup, Method Editor, Maintenance, and Configuration. The 'Home' tab is active. On the left side, there is a vertical menu with options: Overview, Graphs, and Queue. The 'Queue' option is selected and highlighted in blue. The main area of the interface shows a table with the following data:

#	Auto-continue	Submitter	Job name	Duration
	Ready	admin	6 x BSA 30min	11:04:15
1	<input type="checkbox"/>	admin	admin-20100824-1714	11:04:15

Below the table, there is a large empty rectangular area. At the bottom of the interface, there is a toolbar with several buttons: Exit..., Delete, Up, Down, Edit, and Properties. On the left side of the main area, there is a 'Job' label with a value of '00:00:00'. Below that, the date and time '24 aug 2010 17:14' and the user name 'admin' are displayed.

The Job Properties box for the Sample list appears (see [Figure 44](#)).

Figure 44. Sample list job properties on the Queue page

The screenshot shows the Thermo Scientific software interface. At the top, there are navigation tabs: Home, Batch Setup, Method Editor, Maintenance, and Configuration. On the left, there is a sidebar with buttons for Overview, Graphs, and Queue. The main area displays a table of jobs in the queue. A pop-up window titled 'admin-20100825-0002 job properties' is open, showing a table of job details.

#	Auto-continue	Submitter	Job name	Duration
	Ready	admin	6 x BSA 30min	11:04:15
1	<input type="checkbox"/>	admin	admin-20100824-1714	11:04:15

	Position	Name	Method
1	0-A1	Sample 1	BSA 30 min RePlay
2	0-B1	Sample 2	BSA 30 min RePlay
3	0-D1	Sample 4	BSA 30 min RePlay
4	0-E1	Sample 5	BSA 30 min RePlay
5	0-F1	Sample 6	BSA 30 min RePlay

At the bottom of the pop-up window, there are buttons for 'Close' and 'Show batch'. In the background, the Queue page shows a 'Job' with a duration of '00:00:00', a timestamp of '24 aug 2010 17:14', and the user 'admin'. There are also 'Exit...' and 'Delete' buttons at the bottom of the Queue page.

IMPORTANT Reschedule the batch after you complete the batch edit; otherwise, the instrument stops after the current sample run.

Troubleshooting a Sample Run

If a problem occurs during a sample analysis, the EASY-nLC II instrument displays the error in a pop-up box and stops the sample processing.

In this situation, the safest approach is to press Cancel Batch and then go to Maintenance > Scripts (see [“Executing Maintenance Scripts”](#) on page 46). Here you can try to localize/fix the error through maintenance scripts/jobs (see [“Maintenance Scripts”](#) on page 105).

When the problem is fixed, reschedule the batch (possibly changing the number of samples, their well positions, or both).

Autosampler Calibration

This chapter describes how to calibrate the autosampler.

Use this procedure in the following situations:

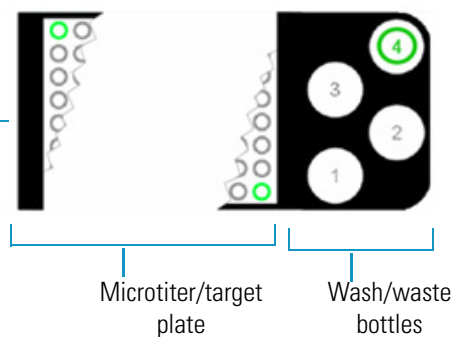
- When you want to use a new type of target plate or Microtiter plates
- When you would like to replace the needle
- For problems with the sample pickup: The needle is not placed in the center of the well, or not all of the liquid is picked up from one or more of the vials.
- For problems with the wash/waste containers: The needle is not placed in the center of the bottle.

Contents

- [Replacing the Autosampler Adapter Plate](#)
- [Managing Plate Formats](#)
- [Preparation and Basic Principles](#)
- [Calibrating Plates](#)
- [Calibrating the Wash/Waste Compartment](#)

The calibration covers the following areas of the autosampler:

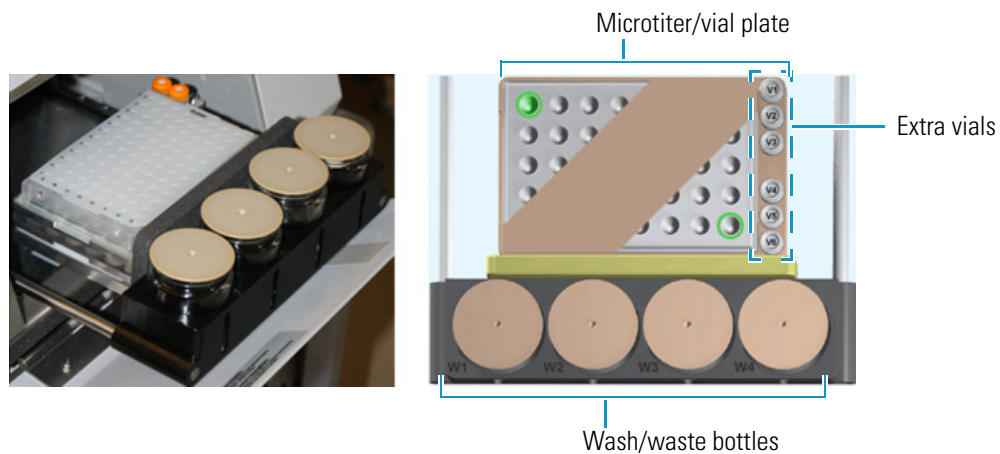
- Original autosampler (ASA)



7 Autosampler Calibration

Replacing the Autosampler Adapter Plate

- Latest autosampler (ASC)



Replacing the Autosampler Adapter Plate

Depending on your autosampler type, follow the appropriate procedure:

- To replace the adapter plate on the ASA autosampler
- To replace the adapter plate on the ASC autosampler

❖ To replace the adapter plate on the ASA autosampler

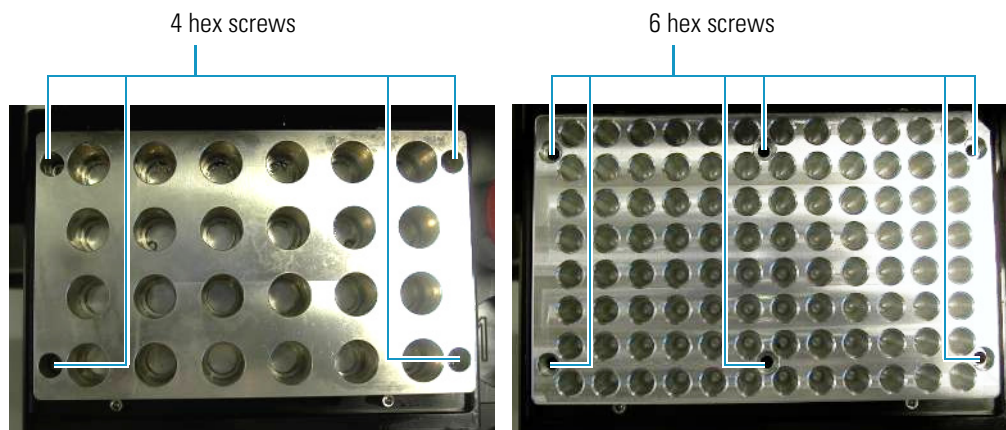
1. Go to **Home > Overview** to eject the tray.



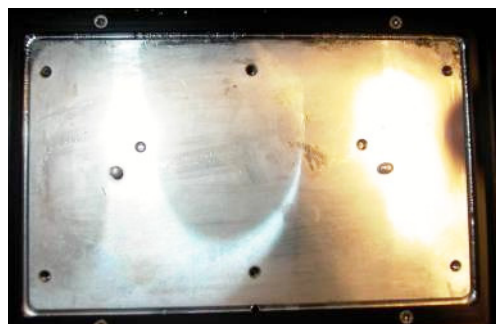
2. Use the 2.5 mm hex key provided with your system to remove the current plate from the autosampler tray. Turn the hex screw counterclockwise.



The adapter plate is mounted with 4 or 6 hex screws:



3. Place the new adapter plate on the tray.



4. Fasten the hex screws in the following balanced order:

- a. Screw in the upper right corner.
- b. Screw in the lower left corner.
- c. Screw in the upper left corner.
- d. Screw in the lower right corner.
- e. Screws in the middle.

Note Mounting the screws in a balanced order can avoid distortion of the adapter plate.

The autosampler is now ready for new plate formats. First, you must calibrate the Microtiter/target plates. See [Managing Plate Formats](#), [Preparation and Basic Principles](#), and [Calibrating the Wash/Waste Compartment](#).

7 Autosampler Calibration

Replacing the Autosampler Adapter Plate

❖ To replace the adapter plate on the ASC autosampler

1. Use the Overview page of the Home menu to eject the tray.



2. Remove the old adapter plate.



3. Place the new adapter plate in the autosampler.

The autosampler is now ready for new plate formats. First, you must calibrate the Microtiter/target plates. See [Managing Plate Formats](#), [Preparation and Basic Principles](#), and [Calibrating the Wash/Waste Compartment](#).

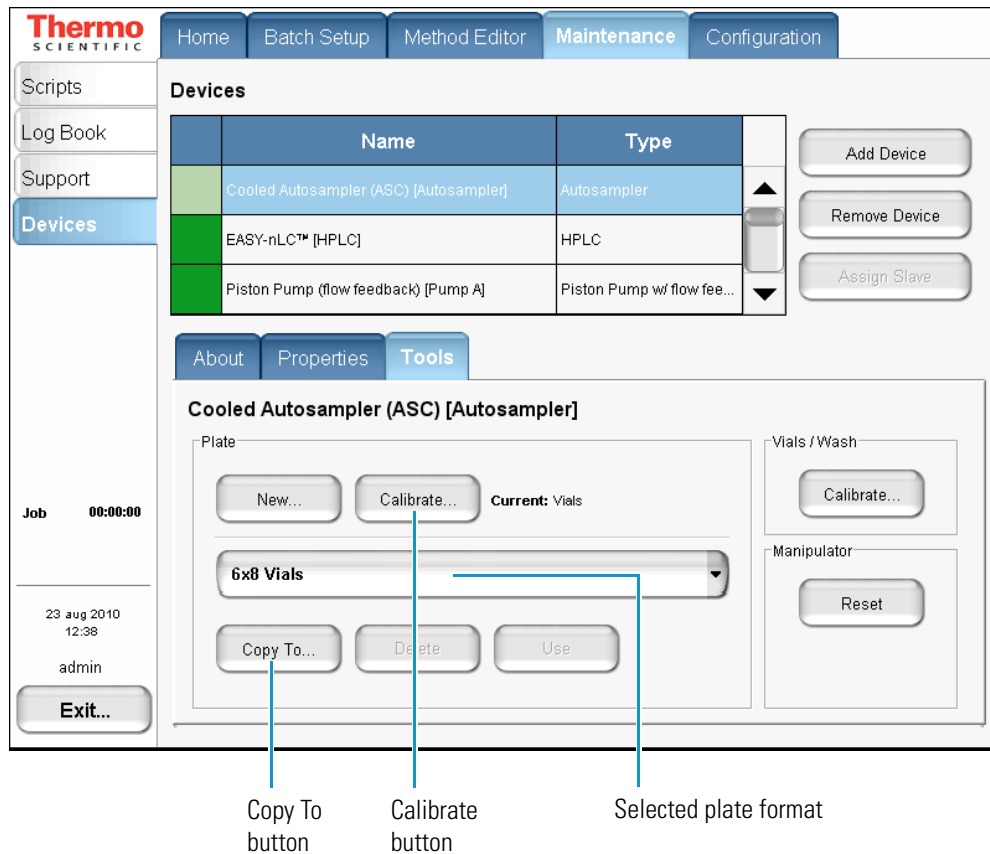
Managing Plate Formats

Follow the appropriate procedure to use an existing plate format, create your own plate format, or delete a format you no longer need.

❖ **To create a new plate format**

1. Go to **Maintenance > Devices**. See [Figure 45](#).

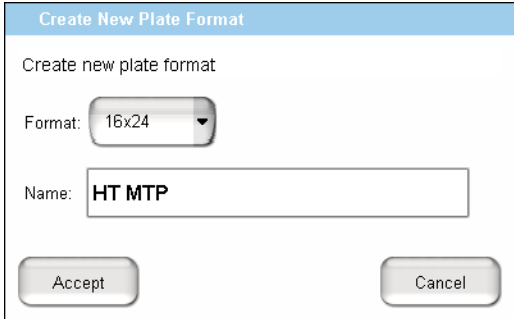
Figure 45. Tools page under Maintenance > Devices showing plate format selection to copy



2. In the Devices list, select **Cooled Autosampler** and then press the **Tools** tab.

3. Do one of the following:
 - a. Press **New** to open the Create New Plate Format dialog box (Figure 46).

Figure 46. Create New Plate Format dialog box



- b. Select a format from the Format list and type in a name.
- c. Press **Accept**.
- d. Under the Tools tab, press **Calibrate**.

–or–

- On the Tools page, select an existing plate format from the Plate list and press **Copy To** to create a new plate format based on your selection (see Figure 45).

❖ **To select a given plate format**

1. Go to **Maintenance > Devices** (see Figure 45).
2. In the Devices list, select **Cooled Autosampler** and then press the **Tools** tab (see Figure 45).
3. Under Tools, select a plate format from the Plate list.
4. Press **Use**.

Note The plate selection in your MS control software must match the instrument selection.

❖ **To delete a plate format**

1. Go to **Maintenance > Devices** (see Figure 45).
2. In the Devices list, select **Cooled Autosampler** and then press the **Tools** tab (Figure 45).
3. Under Tools, select a plate format from the Plate list.
4. Press **Delete**.

Note You cannot delete the current plate.

Preparation and Basic Principles

Explanation of autosampler coordinates:

<i>x</i> axis	On the Microtiter plate: Direction A to H
<i>y</i> axis	On the Microtiter plate: Direction 1 to 12
<i>z</i> axis	Needle height

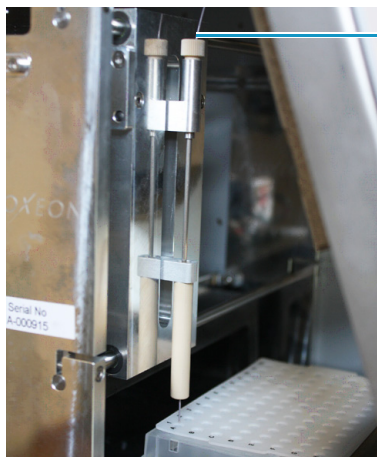


CAUTION Calibration of the autosampler requires removal of the side plate and visual inspection of the autosampler needle holder. Because the autosampler compartment contains moving parts and sharp needles, make sure to keep hands clear when operating the autosampler during calibration.

❖ To calibrate vials and microtiter plates

1. Ensure that the microtiter plates are in place in the tray.
2. Go to **Maintenance > Devices**.
3. From the Devices list, select the autosampler.
4. Press the **Tools** tab below the Devices list.
5. In the Plate area, select a plate format from the list, or create a new plate format.
(To create and select plate formats, see “[Managing Plate Formats](#)” on [page 75](#).)
6. Press **Calibrate** to calibrate the selected plate format. The Calibration dialog box opens.
7. In the Calibration dialog box, press **Eject** to eject the autosampler tray.
8. Check that the correct adapter plate is mounted on the autosampler.
9. Place a Microtiter plate on the adapter plate and cover the Microtiter plate with a plastic film or a rubber mat.
10. Insert the autosampler tray and press **Insert**.

11. Remove the left side panel so that you can watch the movement of the needle.



Use a black marker pen to make a line on top of the PEEKsil tubing.

12. Calibrate the plate. (See “Calibrating Plates,” next.)

Calibrating Plates

For information on selecting an existing plate or creating a new one, see “Managing Plate Formats” on page 75.

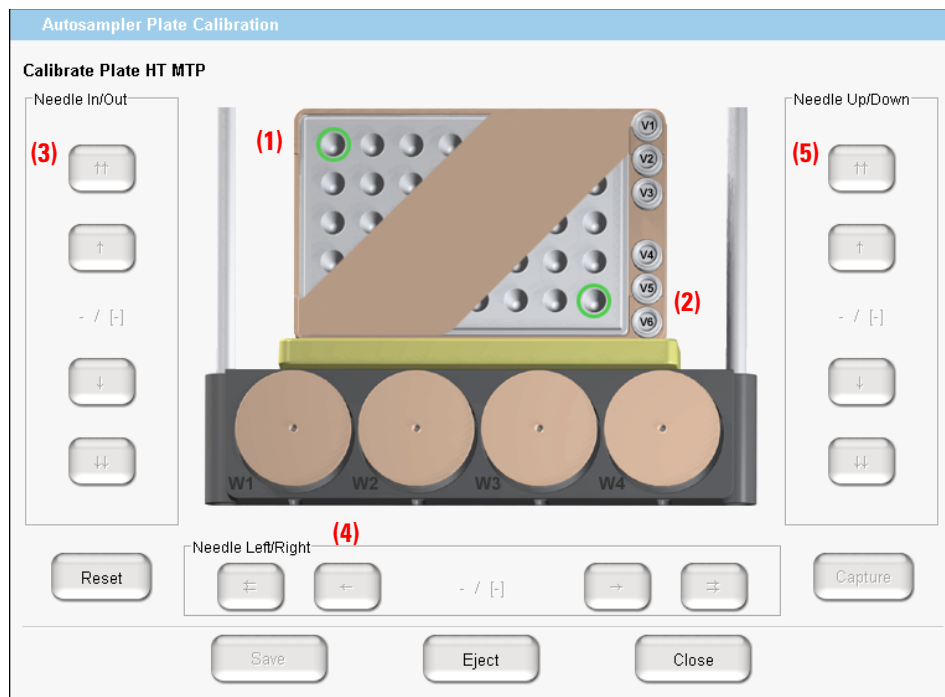
IMPORTANT When calibrating plates on the latest autosampler (ASC), make sure to remove the vial caps or plate covers. Keeping the caps or covers on makes the needle susceptible to bending.

❖ To calibrate plates

1. Put empty vials with no caps in the top left and bottom right positions (position A1 (1) and F8 (2) in the case of the default 6 × 8 vial adapter shown in Figure 47).
2. Go to **Maintenance > Devices**.
3. From the Devices list, select the autosampler.
4. Under the Tools tab in the Plate area, press **Calibrate** (see Figure 45 on page 75). The Autosampler Plate Calibration dialog box opens (Figure 47).

5. Press the green well in the upper left corner of the plate view (1). This moves the needle to the current saved coordinate.

Figure 47. Plate view in the Autosampler Plate Calibration dialog box



Note The EASY-nLC II application copies the saved coordinate to the left side of the slash (/), except for Needle Up/Down, which starts with the top position. The left side coordinate is the working/current coordinate and is not saved yet.

6. Center the needle above the well by pressing the **Needle In/Out** (3) and **Needle Left/Right** (4) buttons.
7. Lower the needle by pressing the **Needle Up/Down** buttons (5) until the needle reaches a level of 2–3 mm (0.08–0.12 in.) above the well. Start with the big step button (⇕) and change to the small step button (↓) when you are near the well.

8. Move slowly with the small step button (↓) and lower the needle until it reaches the bottom of the well.

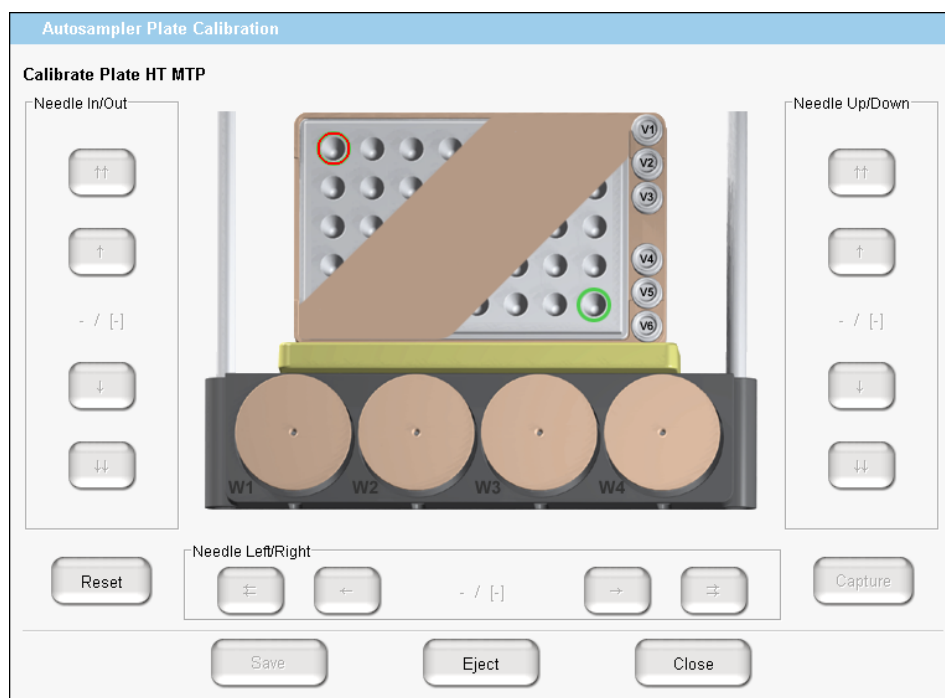


Observe when the black mark on the PEEKsil tubing moves up a bit.

Note The PEEKsil at the top of the needle holder rises a little when the needle reaches the bottom of the well. You can observe this when the black mark moves.

9. When you are satisfied with the current *xyz*-coordinate for the well bottom, press **Capture**. The color of the well changes from green to red (see [Figure 48](#)).

Figure 48. Plate calibration captured



10. Press the green well in the lower right corner of the plate view (see (2) in [Figure 47](#)), which moves the needle to the current saved coordinate.
11. Repeat [step 6](#) to [step 9](#) on that well.

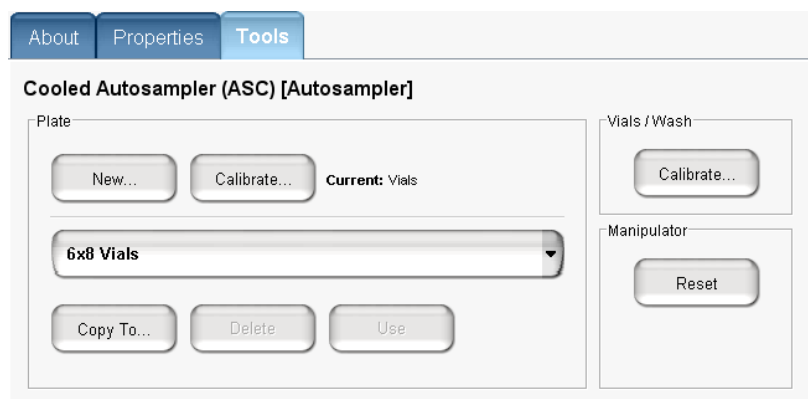
12. Press **Save** and then press **Close** to close the calibration dialog box.
13. Refit the side plate on the system.
14. Replace vial caps and plate covers as necessary.

Calibrating the Wash/Waste Compartment

Make sure that you have placed an empty vial in position V1 and a wash bottle with insert in position W4.

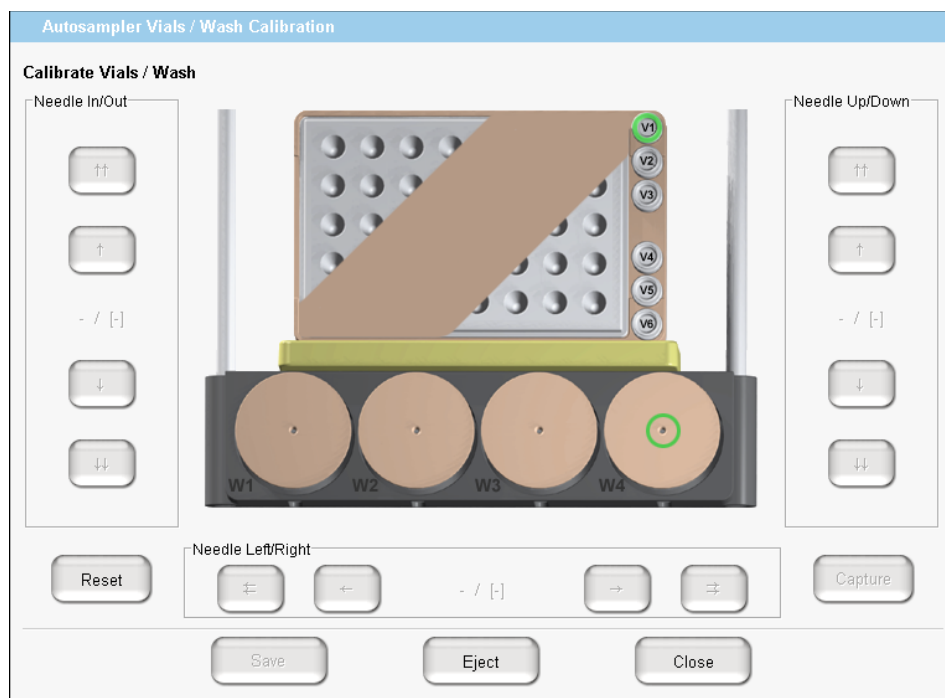
❖ To calibrate the wash/waste compartment

1. Go to **Maintenance > Devices**.
2. From the Devices list, select the autosampler.
3. Press the **Tools** tab below the Devices list.

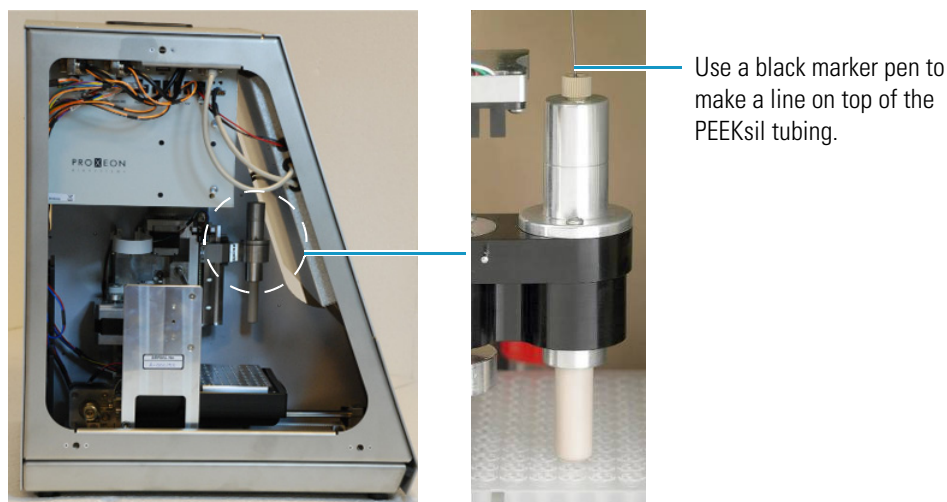


4. In the Vials/Wash area, press **Calibrate**. The Autosampler Vials/Wash Calibration dialog box opens where you calibrate the wash or vial position (see [Figure 49](#) on [page 82](#)).

Figure 49. Autosampler Vials/Wash Calibration dialog box



5. Prepare the bottle for position 4 (used for ejecting waste and cleaning the outside of the injection needle). See “Autosampler Bottles” on page 44.
6. Place the bottle in position W4.
7. Eject the autosampler by pressing **Eject** and place the vial bottle in V1.
8. Insert the autosampler tray if ejected and press **Insert**.
9. Remove the left side panel so that you can watch the movement of the needle.



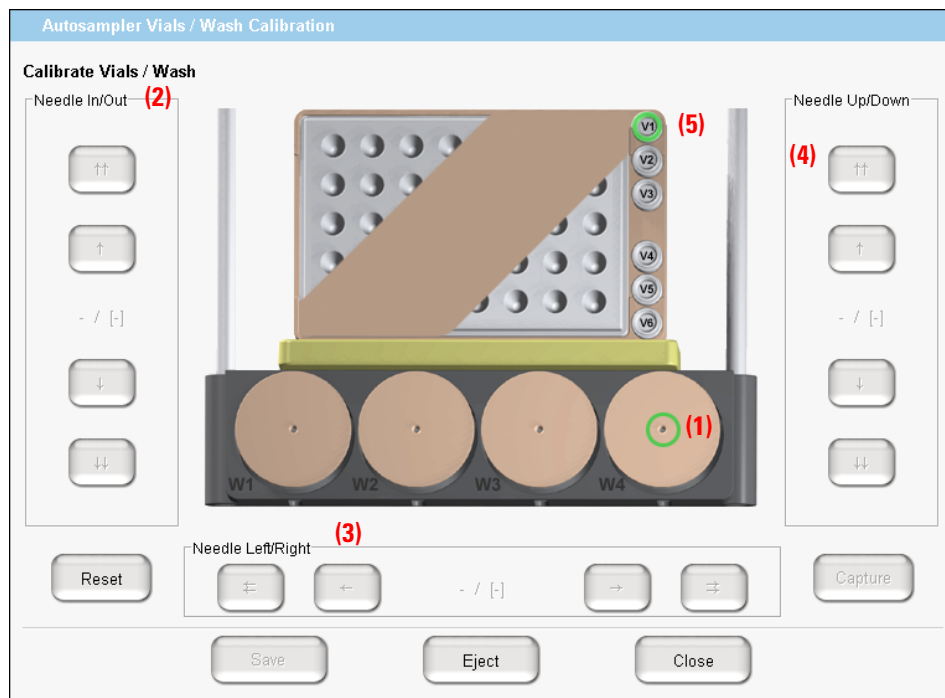
10. Continue with calibrating the wash/waste/vial positions.

❖ **To calibrate the wash/waste/vial positions**

1. Press the bottle container in W4 (1), as show in Figure 50. This moves the needle to the current saved coordinate.

Note The EASY-nLC II application copies the saved coordinate to the left side of the slash (/) except for Needle Up/Down, which starts with the top position. The left side coordinate is the working/current coordinate, which is not yet saved.

Figure 50. Vials/Wash view in the Autosampler Vials/Wash Calibration dialog box



2. Center the needle above the well by pressing the **Needle In/Out** (2) and **Needle Left/Right** (3) buttons.
3. Lower the needle by pressing the **Needle Up/Down** buttons (4) until it reaches a level of 2–3 mm (0.08–0.12 in.) above the bottle. Start with the big step button (⇓) and change to the small step button (↓) when you are close to the well.
4. Using the small step button (↓), slowly let the needle enter the hole and lower the needle until it reaches the bottom of the wash insert in the bottle.

The ASA autosampler has a limit of 850 steps on the z axis (vertical). This limit is set to avoid bent needles. You might be able to go for more than 850 steps during calibration, but the needle does not go more than 850 steps during normal work.

Note The PEEKsil at the top of the needle holder rises a little when the needle reaches the bottom of the well.

7 Autosampler Calibration

Calibrating the Wash/Waste Compartment

5. Press **Capture** when you are satisfied with the current *xyz*-coordinate for position W4. The color of the bottle container number changes from green to red.
6. If the vial needs to be calibrated, press **V1** vial (5) to select it.
7. If needed, repeat [step 2](#) through [step 5](#) for vial A.
8. Press **Save** and then press **Close** to close the calibration dialog box.

❖ To check sample pickup

1. Identify the positions of the wells that you have problems with.
2. Fill the wells with a given volume of solvent A. This example uses 12 µL.
3. Put a plastic film on the Microtiter plate.
4. Insert the tray.
5. On the Maintenance > Scripts page, run the Sample Pickup script with a pickup volume of 10 µL and your default flow (see [Figure 51](#)). Run the script for each well you pointed out in step 1.

Figure 51. Sample Pickup script under Maintenance tab

The screenshot displays three panels of the Sample Pickup script interface. Each panel has a 'Category' dropdown set to 'Test' and a 'Name' dropdown set to 'Sample pickup'. Below these are tabs for 'Description', 'Parameters', and 'Output'.

Panel 1 (Left): The 'Description' tab is active, showing the text: "Test autosampler sample pick up. Please fill plate well with a defined volume before running test. Max volume: Loop volume - 2µl. Max flow: 40.0 µl/min." The 'START' button is highlighted in green.

Panel 2 (Middle): The 'Parameters' tab is active, showing a table with the following data:

Parameter	Value
Volume [µl]	10.00
Flow [µl/min]	20.00
Position	0-A1

The 'START' button is highlighted in green.

Panel 3 (Right): The 'Output' tab is active, showing the text: "'Sample pickup' started Prime pump - filling...". The 'STOP' button is highlighted in red.

6. After the EASY-nLC II application finishes the script:
 - a. Eject the tray.
 - b. Check all wells one by one with a pipette.
 - c. Check that there are 2 μL left in all wells:
 - If there are more than 2 μL left in *all* wells, there might be a problem with the autosampler calibration, or the autosampler tray/adaptor plate is not aligned. Calibrate the autosampler again and do the sample pickup check.
 - If there are more than 2 μL left in some of the wells, there might be air in the system. Run the Flush Air script on pump S (see [Figure 52](#)) and then the sample pickup check.

Figure 52. Flush Air script under Maintenance tab

Category: Prepare Name: Flush air

Description Parameters Output

Flush air from pumps. Make sure to check solvent levels before running.

Parameter	Value
Flush pump A	<input checked="" type="checkbox"/>
Flush pump B	<input checked="" type="checkbox"/>
Flush pump S	<input checked="" type="checkbox"/>
Flush vol. threshold [μL]	10.00

'Flush air' started
Pump A: Iteration 1
Pump B: Iteration 1
Pump B: Refilling...
Pump A: Refilling...

Schedule START STOP

You are finished when there are 2 μL left in all wells.

Troubleshooting

This chapter offers help with the EASY-nLC II system in the form of troubleshooting tips. The following tables provide troubleshooting tips in several different areas.

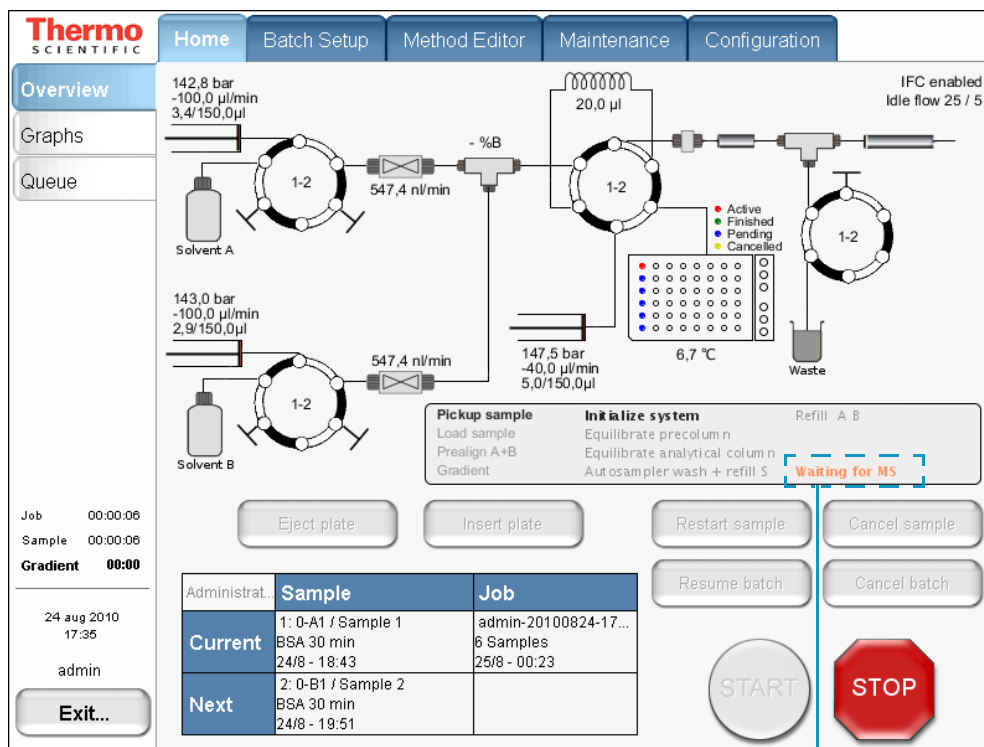
Contents

- [Running a Batch](#)
- [Autosampler](#)
- [Contact Closure](#)
- [Network Access](#)
- [Starting Up the System](#)
- [System Component States \(Valves, Pumps, Autosampler, and Plate Cooler\)](#)
- [Miscellaneous](#)

Running a Batch

For troubleshooting tips when running a batch, see [Table 12](#).

Figure 53. Waiting for MS contact closure signal on the Overview page



Message in activity box

Table 12. Running a batch troubleshooting tips (Sheet 1 of 2)

Symptom	Possible causes	Action
Gradient does not start.	Waiting for contact closure (see Figure 53)	Check the contact closure connections.
Maximum pressure is exceeded during sample loading or reequilibration (IFC disabled), or pressure is maintained but the flow is dramatically decreased (IFC enabled).	Possible blocked column	<ol style="list-style-type: none"> 1. Remove all columns from the system. 2. Run the Precolumn Equilibration script with a flow of 10 $\mu\text{L}/\text{min}$. This should give you a back pressure of approximately 80 bar.

Table 12. Running a batch troubleshooting tips (Sheet 2 of 2)

Symptom	Possible causes	Action
Max pressure is exceeded during gradient.	Back pressure that is too high on the column for the specified flow	Reduce the flow in the method and schedule a new batch with the changed method.
	Blocked flow paths	Contact Thermo Fisher Scientific for advice: easysupport@proxeon.com
It takes extremely long before flow rate stabilizes.	Air in the system	Run the maintenance script: Purge/Flush All See “Purging and Flushing the Pumps” on page 47.
	Leakage in the system	Run the maintenance script: “Test – Leaks” on page 113
XYZ robot error	Possible step loss on one of the motors because of a blockage in the movement of the axis	Remove the obstacle and try again. You might need to calibrate the autosampler for more precise penetration of the Microtiter plate or wash/waste containers.
	Repetition of the error without any visible reason	Contact Thermo Fisher Scientific: easysupport@proxeon.com
Time-out on flow sensor	Air in the pumps	Run the maintenance script: “Prepare – Flush Air” on page 107
Sample is not being picked up.	Air in the pumps	On the ASA: Check that the solvent level in W3 is more than 2/3 full; otherwise, refill it.
		Run the maintenance script: “Prepare – Flush Air” on page 107 If this does not work, try running the Leak Test on pump S (see “Test – Leaks” on page 113) and then the Flush Air script.
Valve – unknown position	Valve trimming needed	Run the maintenance script: “Test – Valve Check” on page 114

Autosampler

See [Table 13](#) for tips on handling autosampler issues. To access the Autosampler control window ([Figure 54](#)), go to **Home > Overview**, and press the plate icon.

Figure 54. Autosampler control window

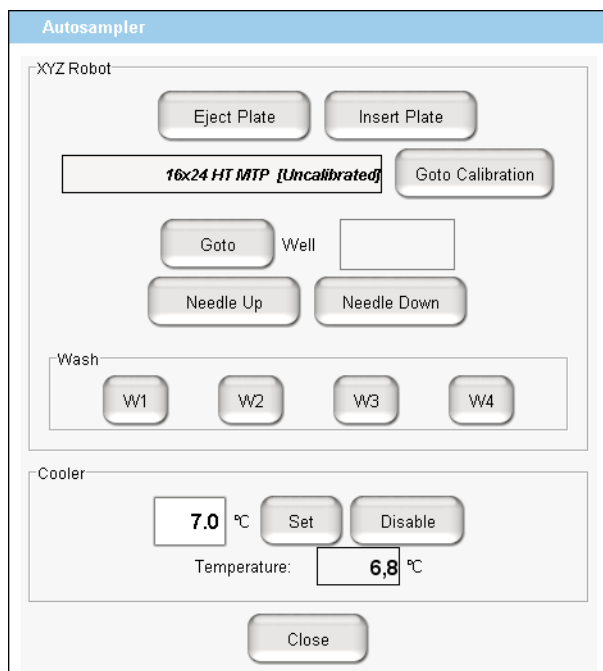
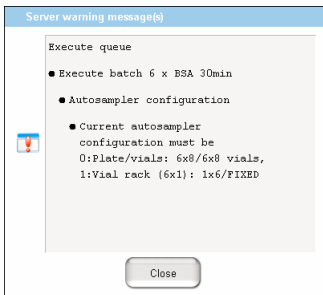


Table 13. Autosampler troubleshooting tips (Sheet 1 of 2)

Symptom	Possible causes	Action
Sample pickup fails.	Bad autosampler calibration or air in the sample (S) pump	<ol style="list-style-type: none"> 1. Fill some of the wells with a given volume of solvent A, for example, 12 μL. 2. Put a plastic film on the Microtiter plate. 3. Run the maintenance script: “Test – Sample Pickup” on page 112, with a pickup volume of 10 μL and your default flow. Run the script for each well. 4. Eject the tray and check the wells one by one with a pipette. 5. Run the “Prepare – Flush Air” on page 107 script, or calibrate the autosampler if the problem still appears.

Table 13. Autosampler troubleshooting tips (Sheet 2 of 2)

Symptom	Possible causes	Action
The needle cannot penetrate the plastic film on the Microtiter plate.	Damaged or bent needle tip	<ol style="list-style-type: none"> 1. Move the autosampler to position A1 by using the Autosampler control window on the Home > Overview page. See Figure 54 on page 90. 2. Remove the left side panel and gently push the PEEKsil cover upwards on the autosampler assembly to expose the PEEKsil needle. The needle should not be bent or damaged in any way. 3. If there is any residue on the needle tip, use tissue with some ethanol to wipe it. 4. Let go of the PEEKsil cover again and check if the PEEKsil needle is on level with the cover (the PEEKsil cover should completely cover the needle).
Cannot run a batch. The following or similar error message pops up when you run the batch:	<p>Plate format deleted from your configuration with the Delete button from Maintenance > Devices > Autosampler > Tools</p> 	If you want to run the batch, you must create the plate format again.
The needle cannot reach the bottom of the solvent bottles (position 1–3).	The needle in the original autosampler (ASA) is not designed to reach the bottom of the solvent bottles.	None.

Contact Closure

Symptom	Possible causes	Action
<p>The MS is waiting for a contact closure signal.</p> <p>-or-</p> <p>The EASY-nLC II system is waiting for a ready signal from the MS.</p>	Wrong contact closure setup	<p>If you are running two-way contact closure (with feedback from the MS), set the contact closure Protocol setting to Two-way on the Configuration > Connections page.</p> <p>If you are running one-way contact closure (that is, the EASY-nLC II system sends a start signal to the MS, but feedback from the MS is ignored), set the contact closure Protocol setting to One-way on the Configuration > Connections page.</p>

Network Access

Symptom	Possible causes	Action
System is not responding after pressing Save Configuration on the Network page under Configuration.	Length of completion time to update the network settings	As indicated, allow for wait time while the system changes network settings.
Cannot access the EASY-nLC II system through the network.	Changed network address	Go to Configuration > Network and press Save Configuration .

Starting Up the System

Symptom	Possible causes	Action
Dark monitor	Broken or open main power fuse	Remove the right side panel of the EASY-nLC II instrument. If there is no light on the pumps, switch off the EASY-nLC II instrument and check the main fuse.
	Defective monitor	Attach an external monitor to the VGA connector at the back of the instruction (labeled MONITOR).
Startup bar is no longer proceeding.	Damaged system	Switch off the EASY-nLC II instrument and wait five seconds before switching it on again. In some cases this recovers the system. Contact Thermo Fisher Scientific if the problem recurs: easysupport@proxeon.com

System Component States (Valves, Pumps, Autosampler, and Plate Cooler)

Symptom	Possible causes	Action
The LED on the device is blinking red.	Device in bootloader mode	The device stopped working. This could also block the function of other devices. Contact Thermo Fisher Scientific: easysupport@proxeon.com

Miscellaneous

For additional issues that might arise while using the EASY-nLC II system, consult these miscellaneous troubleshooting tips (Table 14).

Table 14. Miscellaneous troubleshooting tips (Sheet 1 of 2)

Symptom	Possible causes	Action
Copying log files to a USB memory stick failed.	Memory stick in wrong format	Format the memory stick in FAT/FAT16 format.
	Memory stick not recognized by the system	Use one of the memory sticks supplied by Thermo Fisher Scientific, or try plugging the stick into the other USB port. Contact Thermo Fisher Scientific if neither action works: easysupport@proxeon.com
Copying to the USB memory stick does not stop.	Too much data to be copied	Remove the memory stick from the EASY-nLC II system and press Ignore when the error message appears. Contact Thermo Fisher Scientific for a system check-up: easysupport@proxeon.com
Forgot the admin password.		If you are already logged into the EASY-nLC system, Thermo Fisher Scientific can extract the password from the system in two ways: <ul style="list-style-type: none"> • If the network is set up for remote access, choose Maintenance > Support and press Connect. –or– • Run a factory reset.

Table 14. Miscellaneous troubleshooting tips (Sheet 2 of 2)

Symptom	Possible causes	Action
System is running slower and slower.	Memory filled with graph data from long gradient run	Choose Home > Graph and disable the graphs by pressing No Graph in each graph window.
	System overload caused by other reasons	Restart the application or power down/power up from Maintenance > Scheduling .
Unhandled error or SVG (scalable vector graphics) error	Program error in the software release	<p>In most situations you can continue your work by pressing Ignore or OK. More serious error situations might require restarting the HPLC.</p> <p>To improve on the software quality, e-mail details to: easysupport@proxeon.com</p> <p>Or, connect to the remote support server. See “Connecting the EASY-nLC II System to the Support Server” on page 97.</p>

Remote Support

The EASY-nLC II system is designed with an integrated remote support function. While connected, Technical Support can follow the processes on your system to better understand any issues that arise.

Contents

- [Connecting the EASY-nLC II System to the Support Server](#)
- [Saving System Files on a USB Removable Storage Device](#)

Connecting the EASY-nLC II System to the Support Server

Check that your EASY-nLC II system is connected to the network before you continue. For more information, see “[Network Connection](#)” on [page 36](#).

❖ To connect to the support server

1. Choose **Maintenance > Support** (see [Figure 55](#)).

The options for the remote connection are in the Connect to Proxeon Support area.

9 Remote Support

Connecting the EASY-nLC II System to the Support Server

Figure 55. Support page under the Maintenance tab

The screenshot shows the Thermo Scientific software interface. At the top, there are navigation tabs: Home, Batch Setup, Method Editor, Maintenance (selected), and Configuration. On the left side, there is a sidebar menu with options: Scripts, Log Book, Support (selected), and Devices. The main content area is titled "Connect to EASY-nLC support:". It includes a "Remote Port Number: 2000" field, a "Connect" button, a "Disconnect" button, and a "Status: Not Connected" box. Below this is a "Message window (can be used to chat with support when connected):" with "Show" and "Clear" buttons. The next section is "Log file copy:", featuring radio buttons for "To home directory" (selected) and "To USB removable storage", a "Copy Log Files" button, and a "Status: Idle" box. The final section is "Contact:", which lists contact information for Thermo Fisher Scientific: Edisonsvej 4, DK-5000 Odense C, Denmark. It also provides the email address easysupport@proxeon.com, the website http://www.proxeon.com/hplc, the phone number +45 6557 2300, and the fax number +45 6557 2301. In the bottom left corner, there is a "Job" section showing "00:00:00", a date and time "30 Aug 2010 14:14", the user "admin", and an "Exit..." button.

2. Press **Connect**. A pop-up box opens.

The pop-up box is titled "Message to support team". It contains the following fields: "Name:" with a text input box, "Email address:" with a text input box, "Phone number:" with a text input box, and "Enter message:" with a larger text area. At the bottom of the box, there are two buttons: "Accept" and "Cancel".

3. In the pop-up box, briefly describe the problem that you are experiencing. Provide your contact information and press **Accept**.

The Status box on the right side of the Support page changes from Not Connected to Connected.

The screenshot shows the "Status:" label next to a text box containing the text "Connected!".

The EASY-nLC II system is now connected to the support server. It stays connected until you press Disconnect or the network connection is broken.

Inform Thermo Fisher Scientific if you would like to initiate remote support, and the Technical Support team will start to monitor and diagnose your system.

Saving System Files on a USB Removable Storage Device

The EASY-nLC II application is designed with network access in mind. Through the network you can back up your system, export and import batches and methods, analyze your system by examining the log files, and perform many other tasks. However, when you are not connected to the network or the network is broken, or you do not have a computer with network access that is near the EASY-nLC II system, you can use the copy function. You can copy all the system files on the system to a USB removable storage device, such as a USB flash drive or memory stick.

Before continuing, you need a USB storage device formatted with the file system FAT/FAT16 and at least 128 MB of free space. The EASY-nLC II system is shipped with a USB flash drive for your convenience.

❖ To copy system files to a USB storage device

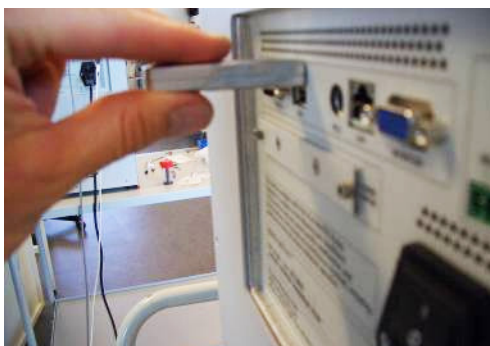
1. Choose **Maintenance > Support**.

The screenshot shows the Thermo Scientific software interface. The top navigation bar includes Home, Batch Setup, Method Editor, Maintenance, and Configuration. The left sidebar contains Scripts, Log Book, Support (highlighted), and Devices. The main content area is titled 'Connect to EASY-nLC support:' and includes a 'Remote Port Number' field set to 2000, 'Connect' and 'Disconnect' buttons, and a 'Status' field showing 'Not Connected'. Below this is a 'Message window' section with 'Show' and 'Clear' buttons. The 'Log file copy:' section has two radio buttons: 'To home directory' (selected) and 'To USB removable storage'. A 'Copy Log Files' button and a 'Status' field showing 'Idle' are also present. The 'Contact:' section provides Thermo Fisher Scientific contact information. At the bottom left, there is a 'Job' status area showing '00:00:00', a date and time '30 Aug 2010 14:14', the user 'admin', and an 'Exit...' button.

9 Remote Support

Saving System Files on a USB Removable Storage Device

2. Insert the USB storage device into the connector panel on the back of the EASY-nLC II instrument. Choose between one of two USB ports on the upper left side of the panel.



3. In the Log File Copy area, select the **To USB Removable Storage** option.
4. Press **Copy Log Files**.

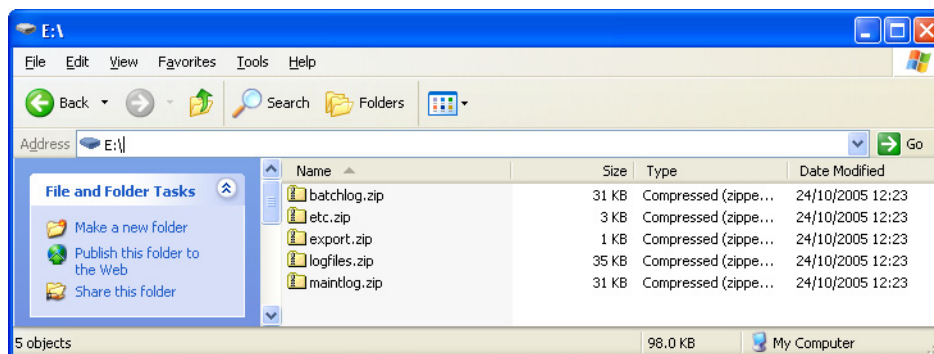
When the copying is finished, the following confirmation appears:



5. Press **Close** and remove the USB storage device from the connector panel.
6. Insert the USB storage device into your computer.

Five zipped files appear on the USB storage device (see [Figure 56](#)).

Figure 56. Windows XP view of the USB memory stick



The files are as follows:

batchlog.zip	Contains one folder for each batch you have run with information on samples, methods, and EASY-nLC II device performance.
etc.zip	Contains system configuration files.
export.zip	Contains exported batches and methods per user and system backup.
logfiles.zip	Contains system log files with information on overall system performance.
maintlog.zip	Contains one folder for each maintenance script you have run with information on EASY-nLC II device performance.

If you have problems with your system, forward the files to Thermo Fisher Scientific Technical Support for assistance.

Maintenance

This chapter provides various maintenance schedules for the EASY-nLC II system, and describes the system's built-in maintenance scripts and when to use them. Consistent use of these procedures will help ensure the longevity of your instrument.

Contents

- [Daily Maintenance](#)
- [Quarterly Maintenance](#)
- [Yearly Maintenance](#)
- [Maintenance Scripts](#)

Daily Maintenance

Check the solvent levels on a daily basis.

❖ To check solvent levels

1. Visually inspect solvent bottle A, solvent bottle B, and the autosampler bottle in position W3 (and the bottles in position W1 and W2 if used). Refill if necessary.
(ASA autosampler only) Make sure that the solvent level in the W1, W2, and W3 autosampler bottles never falls under 2/3 of the bottle volume.

Tip You can also inspect the purity of the solvent to ensure no visible precipitates have formed that might lead to blockages in the pump line.

2. Visually inspect the autosampler wash bottle in position W4 and the waste container in the pump compartment, and empty if necessary.
3. Run the **Purge Solvent** and **Flush Air** scripts with two purge iterations or until flush volumes fall below 10 µL.

Quarterly Maintenance

Perform the following procedures quarterly (every three months).

❖ **To check instrument back pressure**

Run the **Back Pressure** script for both solvents.

❖ **To check the cooler temperature**

Check that the plate temperature is at its set point or approximately 18 to 20 °C (64 to 68 °F) below ambient temperature (readout on the Home > Overview page).

❖ **To check autosampler pickup**

Run the **Sample Pickup** script to check the autosampler pickup function.

❖ **To check the EASY-nLC II system for leaks**

Run the **Leaks** script for “A+B” and for “System,” to check for and pinpoint instrument leaks.

❖ **To check the check valves**

Run the **Purge Solvent** script while checking that solvents are exiting the A and B waste tubing and passing into the waste container when the pumps are ejecting. Make sure that no solvents are going into the pump through these tubes during refill.

Yearly Maintenance

Schedule the following procedures once a year.

❖ **To exchange low pressure solvent filters and high pressure inline filters**

Thoroughly clean new solvent filters before use as these can be a contamination source.

❖ **To check flow sensor calibration**

Run the **Flow Sensor** calibration script with the **Inspection Only** parameter enabled. This checks that the flow sensor calibration is working. Enter **30** minutes for Period.

IMPORTANT Check the valve rotor shifts regularly. When the valve rotor shifts exceed 6000, Thermo Fisher Scientific recommends checking the system for leaks. From this point forward, perform leak tests every 1000 shifts and exchange the rotor when the leak test fails.

Maintenance Scripts

❖ To execute all listed maintenance scripts

1. In the EASY-nLC II application, choose **Maintenance > Scripts**.
2. Select a category and then a name for the specific operation.

You can also schedule some of the scripts for execution using the job queue.

Each of the following sections combines the category with the actual script name.

- Prepare – Purge Solvent
- Prepare – Flush Air
- Prepare – Precolumn Equilibration
- Prepare – Analytical Col Equilibration
- Prepare – Isocratic Flow
- Test – MS Connection
- Test – Sample Pickup
- Test – Leaks
- Test – Valve Check
- Test – Back Pressure
- Test – Autosampler Torque
- Calibrate – Valve Tune
- Calibrate – Flow Sensors
- Calibrate – Reset Pressure Sensor

Prepare – Purge Solvent

The Purge Solvent script fills the chosen pump or pumps with solvent and then ejects it into the waste container. Use the Purge Solvent script when exchanging solvents, removing air from solvent lines, or filling pumps.

❖ To set the Purge Solvent script parameters

1. Press **Purge Iterations** to specify the number of empty/fill cycles you want performed.
2. Select the check boxes for the appropriate pump or pumps.

Note Entering a value of zero (0) for purge iterations will just refill the pumps.

The screenshot displays the 'Maintenance' configuration window for the 'Purge solvent' script. The 'Parameters' tab is active, showing a table with the following data:

Parameter	Value
Purge iterations	10
Purge pump A	<input checked="" type="checkbox"/>
Purge pump B	<input checked="" type="checkbox"/>
Purge pump S	<input checked="" type="checkbox"/>

Below the table are 'START' and 'STOP' buttons, with 'START' being a prominent green button. The right side of the interface features two empty graphs for 'Pump A' and 'Pump B', each plotting 'Flow (nl/min)' and 'Pressure (bar)' against 'Time (min)'. The 'Job' timer shows '00:00:07' and the user is logged in as 'admin'.

Prepare – Flush Air

The Flush Air script pressurizes the pump and then releases that pressure into the flow lines toward the waste container, which is effective for removing air from inside the pump head.

Pressurization time is calculated dynamically while the script is executed and is based on measurements from previous iterations.

The pump pressurizes, releases pressure, and empties for each iteration of the script.

The pressurization is limited to 200 bar and the script measures the amount of pumping required to reach 200 bar. Because the solvents are slightly compressible and the pump itself expands, even an air-free pump allows some pumping before reaching 200 bar. Depending on the solvents, a pumped volume less than 10 μL is acceptable.

❖ To set the Flush Air script parameters

Press **Flush Vol. Threshold** to enter the tolerated threshold for an accepted test result. Below 10 μL is satisfactory.

The script terminates after the first iteration that gives a flush volume below the set threshold.

The screenshot shows the 'Maintenance' configuration window for the 'Flush Air' script. The 'Parameters' tab is selected, showing the following table:

Parameter	Value
Flush pump A	<input checked="" type="checkbox"/>
Flush pump B	<input checked="" type="checkbox"/>
Flush pump S	<input checked="" type="checkbox"/>
Flush vol. threshold [μl]	10.00

At the bottom of the window, there are buttons for 'Schedule', 'START' (a large green button), and 'STOP' (a grey octagonal button). The 'Exit...' button is located in the bottom left corner of the sidebar area.

Prepare – Precolumn Equilibration

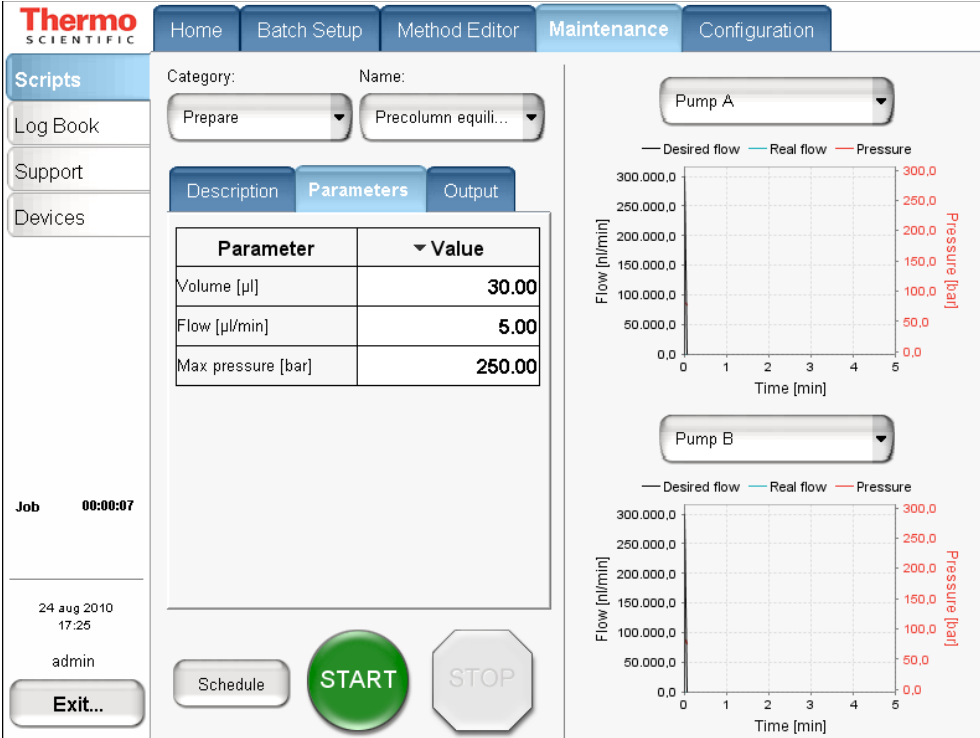
The Precolumn Equilibration script equilibrates the pre-column with solvent from pump A. The script utilizes the IFC system (unless it is disabled), and the procedure below is based on IFC being active. If IFC is disabled, the EASY-nLC II application ignores the Max Pressure setting and the pump is run passively at the specified flow.

Use this script to equilibrate/flush the pre-column, determine a suitable pre-column reequilibration and load flow for method setup, or when initializing a new pre-column.

❖ To set the Precolumn Equilibration script parameters

1. Press **Volume** to set how much solvent A to use for reequilibration of the pre-column. For best results, use 10 column volumes for reequilibration.
2. Press **Flow** to set the flow rate for performing the reequilibration of the pre-column. If the flow field is left empty, the pump operates at the set pressure.
3. Press **Max Pressure** to set the maximum allowed pressure. If the pressure field is left empty, the pump operates at the set flow (as long as it is below the instrument maximum pressure of 300 bar).

If both a flow and a max pressure are specified, the pump flow is limited by whichever parameter is reached first.



The screenshot displays the 'Maintenance' configuration window for the 'Prepare' script. The 'Parameters' section is as follows:

Parameter	Value
Volume [µl]	30.00
Flow [µl/min]	5.00
Max pressure [bar]	250.00

Below the parameters are 'START' and 'STOP' buttons. To the right, there are two empty graphs for 'Pump A' and 'Pump B'. Each graph plots 'Flow (nl/min)' on the left y-axis (0.0 to 300,000.0) and 'Pressure (bar)' on the right y-axis (0.0 to 300.0) against 'Time (min)' on the x-axis (0 to 5). The legend for both graphs indicates 'Desired flow' (black line), 'Real flow' (blue line), and 'Pressure' (red line).

Prepare – Analytical Col Equilibration

The Analytical Col Equilibration script equilibrates the analytical column with solvent from pump A. The script utilizes the IFC system (unless it is disabled), and the procedure below is based on IFC being active. If IFC is disabled, the EASY-nLC II application ignores the Max Pressure setting and the pump runs passively at the specified flow.

Use this script to equilibrate/flush the analytical column, determine suitable analytical reequilibration flow for method setup, or initialize a new analytical column.

❖ To set the Analytical Col Equilibration script parameters

1. Press **Volume** to set how much solvent A to use for reequilibration of the analytical column. For best results, use 10 column volumes for reequilibration.
2. Press **Flow** to set the flow rate for performing the reequilibration of the analytical column. If the flow field is left empty, the pump operates at the set pressure.
3. Press **Max Pressure** to set the maximum allowed pressure. If the pressure field is left empty, the pump operates at the set flow (as long as it is below the instrument maximum pressure of 300 bar).

If both a flow and a maximum pressure are specified, the pump flow is limited by whichever parameter is reached first.

The screenshot displays the Thermo Scientific software interface for the Analytical Col Equilibration script. The interface is divided into several sections:

- Navigation:** Home, Batch Setup, Method Editor, Maintenance (selected), Configuration.
- Scripts:** Log Book, Support, Devices.
- Script Selection:** Category: Prepare, Name: Analytical col eq...
- Parameters Table:**

Parameter	Value
Volume [µl]	10.00
Flow [µl/min]	5.00
Max pressure [bar]	250.00
- Graph:** A line graph showing Flow [nl/min] (left y-axis, 0.0 to 300.000.0) and Pressure [bar] (right y-axis, 0.0 to 300.0) over Time [min] (x-axis, 0 to 5). The graph shows a blue line for 'Desired flow' and a red line for 'Pressure'. The 'Real flow' is shown as a green line. The graph is currently set to 'No graph'.
- Job Information:** Job 00:01:02, 24 aug 2010 23:55, admin.
- Controls:** Exit..., Schedule, START (green button), STOP (grey button).

Prepare – Isocratic Flow

The Isocratic Flow script runs solvent A and B at a fixed mixture ratio. The script utilizes the AFC system to accurately control the flow.

Use this script to tune the mass spectrometer at a given B percentage or when cleaning the instrument/columns.

❖ To set the Isocratic Flow script parameters

1. Press **Volume** to specify the total volume to be delivered by the pumps.
2. Press **Flow** to specify the flow rate to be used.
3. Press **AB Mix** to specify the solvent composition.

The screenshot displays the Thermo Scientific software interface for the 'Prepare' script. The 'Parameters' tab is active, showing a table with columns 'Parameter' and 'Value'. The parameters are: Volume [µl] = 100.00, Flow [µl/min] = 0.30, AB Mix [%B] = 30, and Run indefinitely (checked). To the right, there are two graphs for Pump A and Pump B, each showing 'Desired flow', 'Real flow', and 'Pressure' over a 5-minute period. The 'START' button is highlighted in green.

Parameter	Value
Volume [µl]	100.00
Flow [µl/min]	0.30
AB Mix [%B]	30
Run indefinitely	<input checked="" type="checkbox"/>

Test – MS Connection

The MS Connection script tests the contact closure function between the EASY-nLC II system and the mass spectrometer.

Use this script when testing contact closure.

There are no parameters associated with the MS Connection script. Instead, it uses the settings from the Configuration > Connection page (see “[Setting Up the Mass Spectrometer Connection](#)” on page 35).

For a description of the MS Connection script, see [Figure 57](#).

Figure 57. MS Connection script description



Test – Sample Pickup

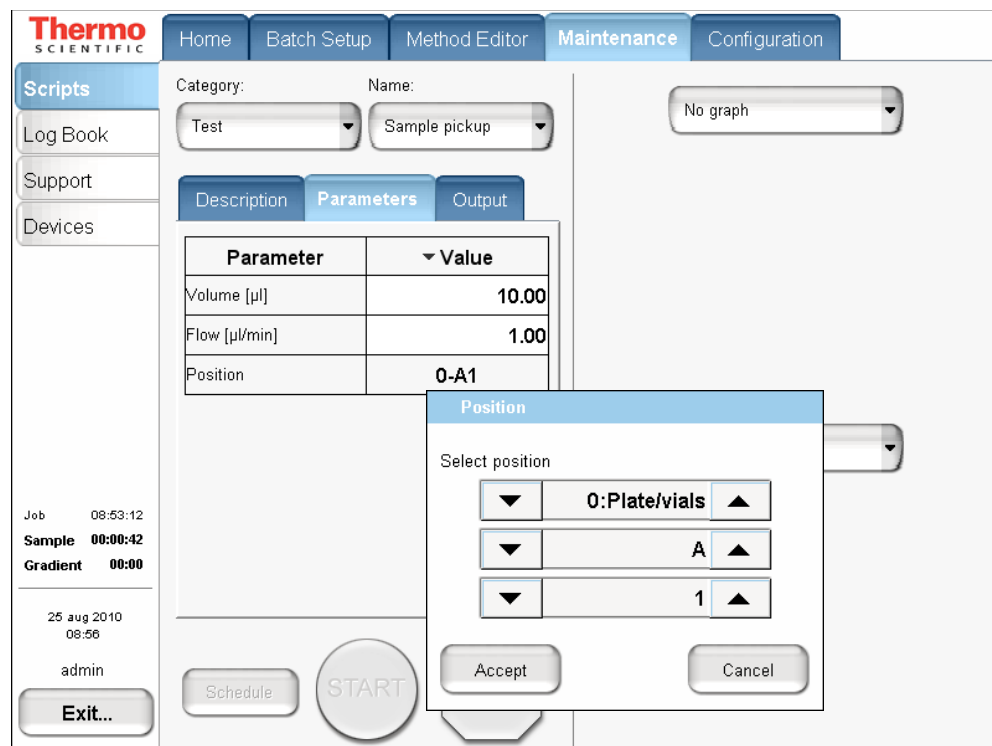
The Sample Pickup script tests the autosampler pickup function by picking up a set volume from the well plate and ejecting it into waste (see [Figure 58](#)).

Use this script when testing the sample pickup function and determining sample pickup flow rates for method setup.

❖ To set the Sample Pickup script parameters

1. Press **Volume** to specify the volume to be picked up.
2. Press **Flow** to specify the flow rate used while aspirating the sample.
3. Press **Position** to specify the position of the sample in the autosampler. See the Position pop-up box in [Figure 58](#).

Figure 58. Sample Pickup script parameters



Test – Leaks

The Leaks script tests for possible leaks on each pump/valve and throughout the flow path, and can also help determine where a possible leak is located. The flow sensors measure the leaks, or they are calculated from pump piston movement.

Use this script when you suspect leak problems in the pumps, the system (flow path), or both.

❖ To set the Leaks script parameters

1. Press the **Parameters** tab.
2. To select which pumps (or system) to perform the leak test on, press the cell in the Value column next to Subsystem(s) (see [Figure 59](#)).

Note You cannot perform the leak test for pump S as it does not operate under pressure during normal batch execution.

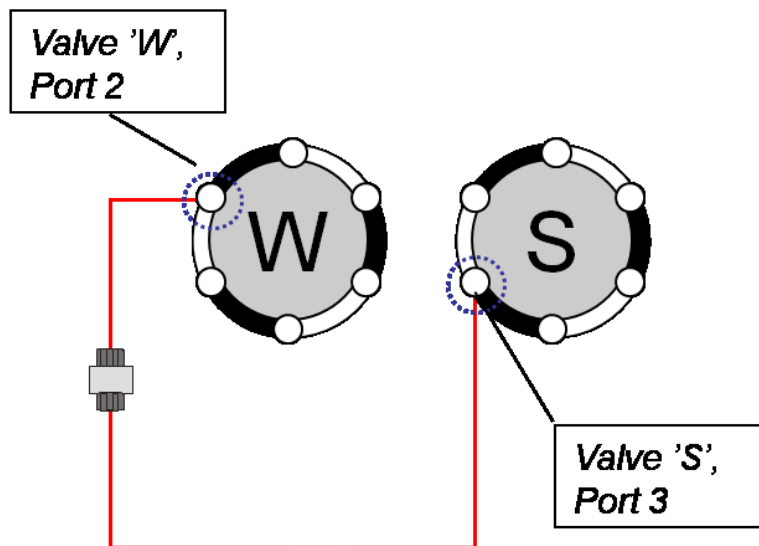
Figure 59. Leaks script parameters

The screenshot displays the Thermo Scientific software interface for configuring the Leaks script. The top navigation bar includes 'Home', 'Batch Setup', 'Method Editor', 'Maintenance', and 'Configuration'. The left sidebar contains 'Scripts', 'Log Book', 'Support', and 'Devices'. The main window is divided into several sections:

- Category and Name:** 'Test' and 'Leaks' are selected in dropdown menus.
- Parameters Tab:** A table with columns 'Parameter' and 'Value'. The 'Subsystem(s)' parameter is set to 'System'. A dropdown menu is open showing options: 'A', 'B', 'A+B', and 'System'.
- Graph:** A line graph showing 'Desired flow' (ml/min) on the left y-axis and 'Pressure' (bar) on the right y-axis against 'Time [min]' on the x-axis. The flow starts at approximately 200 ml/min and drops to 0 at 1 minute. The pressure remains constant at approximately 200 bar. A legend indicates 'Desired flow' (black line) and 'Pressure' (red line).
- Job Information:** 'Job 00:01:02', '24 aug 2010 17:48', and 'admin' are displayed.
- Buttons:** 'Exit...', 'Schedule', 'START' (highlighted in green), and 'STOP' are visible at the bottom.

3. If selecting to perform a System test, ensure that the sample column output line (Valve S, port 3) is connected to waste input (Valve W, port 2) (Figure 60).

Figure 60. Column output line to waste input



Test – Valve Check

The Valve Check script performs a self-test for the selected valves.

Use this script when valves begin to move slowly or cannot reach the required positions and after exchanging valve parts (rotor, stator).

IMPORTANT Only use this script when instructed to by Thermo Fisher Scientific. Make sure to record any numbers that the script generates.

❖ **To set the Valve Check script parameters**

Select the check boxes for the valves that you want to test (Figure 61).

Figure 61. Valve Check script parameters

The screenshot shows the Thermo Scientific software interface for setting script parameters. The 'Maintenance' tab is selected, and the 'Valve Check' script is chosen. The 'Parameters' tab is active, showing a table with the following data:

Parameter	Value
Include valve A	<input checked="" type="checkbox"/>
Include valve B	<input checked="" type="checkbox"/>
Include valve S	<input type="checkbox"/>
Include valve W	<input type="checkbox"/>

Below the table are buttons for 'Schedule', 'START', and 'STOP'. To the right, there are two graphs for 'Pump A' and 'Pump B'. Each graph plots 'Flow [nl/min]' (left y-axis, 0.0 to 300.000.0) and 'Pressure [bar]' (right y-axis, 0.0 to 300.0) against 'Time [min]' (x-axis, 0 to 5). The graphs show 'Desired flow' (black line), 'Real flow' (blue line), and 'Pressure' (red line).

Test – Back Pressure

The Back Pressure script determines instrument back pressure for solvent A and B. The script runs at a preset flow and measures the back pressure on the system.

Use this script when testing for instrument flow path blockage. Before running the script, connect column output to waste input as shown in [Figure 60](#) on [page 114](#).

❖ To set the Back Pressure script parameters

Test solvent A, solvent B, or both A and B by selecting the corresponding check box or check boxes ([Figure 62](#)).

Figure 62. Back Pressure script parameters

The screenshot displays the Thermo Scientific software interface for the Back Pressure script. The interface includes a navigation menu on the left with options like Scripts, Log Book, Support, and Devices. The main area is divided into sections for script configuration and real-time monitoring.

Script Configuration:

- Category: Test
- Name: Back pressure
- Parameters table:

Parameter	Value
Test solvent A	<input checked="" type="checkbox"/>
Test solvent B	<input checked="" type="checkbox"/>

Real-time Monitoring:

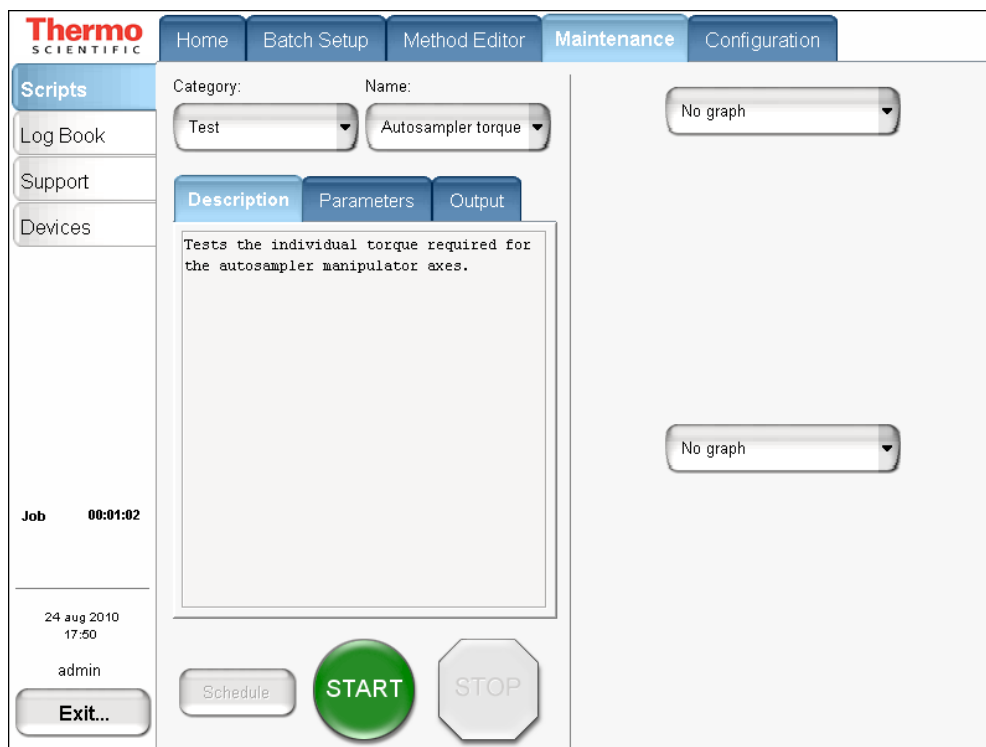
Two graphs are shown, one for Pump A and one for Pump B. Each graph plots Flow (nl/min) on the left y-axis (ranging from -100,000.0 to 300,000.0) and Pressure (bar) on the right y-axis (ranging from 0.0 to 300.0) against Time (min) on the x-axis (ranging from 0 to 5). The graphs show three data series: Desired flow (black line), Real flow (blue line), and Pressure (red line). In both graphs, the desired flow is approximately 200,000.0 nl/min, and the real flow is slightly lower, around 180,000.0 nl/min. The pressure is consistently around 200.0 bar.

Test – Autosampler Torque

The Autosampler Torque script measures the torque required to move the autosampler on each of its axes (Figure 63).

Contact your local Thermo Fisher Scientific representative before running this script.

Figure 63. Autosampler Torque script description



Calibrate – Valve Tune

The Valve Tune script automatically retunes the valve. Do not run this script unless instructed to by a Thermo Fisher Scientific field service engineer.

Calibrate – Flow Sensors

The Flow Sensors script calibrates the flow sensors on pumps A and B.

Use this script when you change the type of solvent in A or B, or you suspect the calibration to be wrong.

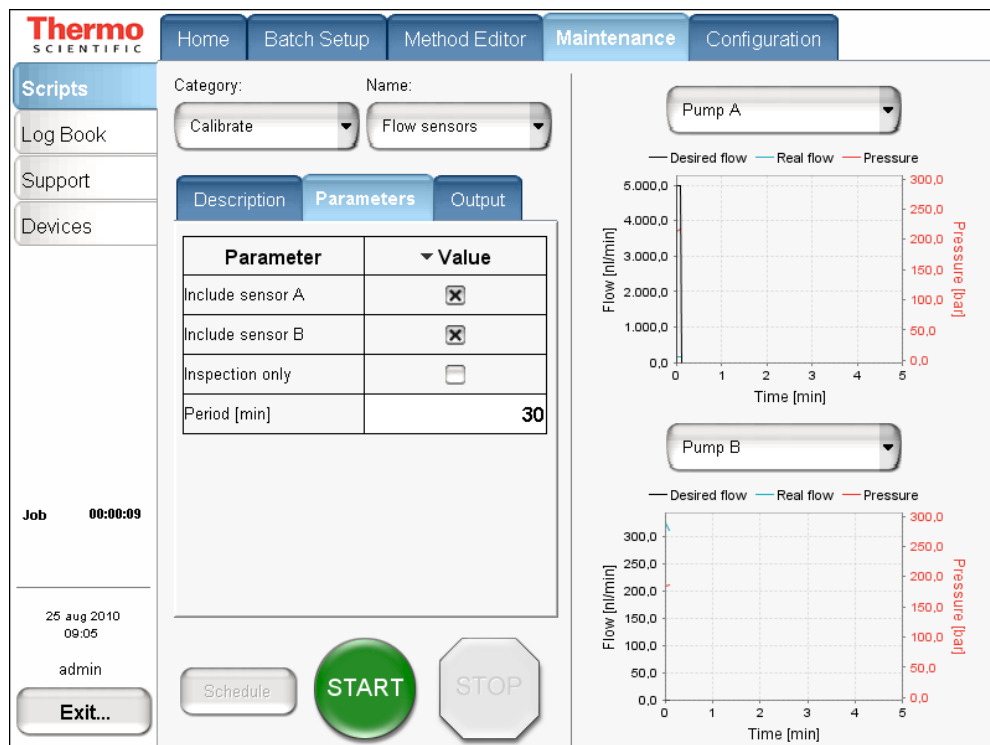
Before calibration, connect a 10 µm ID tubing to the Column Out line to achieve back pressure at approximately 50 bar at 500 nL/min.

❖ **To set the Flow Sensors script parameters**

1. Select the check boxes for the flow sensors that you would like to calibrate (Figure 64).
2. Press **Period** to specify the amount of time (in minutes) for stabilization of the flow before the calibration is performed.

The script makes a 2-point calibration and uses the stabilization period at both calibration points.

Figure 64. Flow Sensors script parameters



Calibrate – Reset Pressure Sensor

This script auto-zeroes the pressure sensor. Contact your local Thermo Fisher Scientific Technical Support representative before running this script.

Specifications

This appendix contains the following specifications for the EASY-nLC II system as well as a flow line schematic.

Contents

- [Performance Specifications](#)
- [Technical Specifications](#)
- [Flow Line Diagram](#)

Performance Specifications

See [Table 15](#) for performance specifications.

Table 15. Performance specification (Sheet 1 of 2)

Item	Specification
Flow range (gradient)	20–2000 nL/min Recommended: 100–1000 nL/min
Flow while loading and equilibrating	Up to 100 μ L/min
Pressure range	0–300 bar
Retention time reproducibility (RT RSD)	Recommended flow range: 0.1 to 0.4% Typically better than 1% outside recommended flow range
Peak widths	Typically 3–5 s fwhm, using EASY-column™ nano-bore columns as supplied at delivery
For more information, go to www.proxeon.com/productrange/nano_lc_easy_columns/	
Carryover	Typically < 0.05% Conditions: Injection of 100 fmol tryptic digests of BSA on EASY-columns as supplied at delivery and by using the Standard Wash procedure for the autosampler
Autosampler pickup volume range	100–18000 nL, with standard loop 100–58000 nL, with up to 60 μ L loop

Table 15. Performance specification (Sheet 2 of 2)

Item	Specification		
Injection reproducibility (injection RSD)	0.2% at 5 μ L pickup 3.0% at 100 nL pickup		
Injection linearity	0.9985 at 0.5–10 μ L injection volume		
Autosampler formats	Original autosampler (ASA):	Latest autosampler (ASC):	
	4 \times 6 Vials	6 \times 8 Vials	+ 6 Vials
	1 \times 96 well MTP	1 \times 96 well MTP	+ 6 Vials
	1 \times 384 well MTP	1 \times 384 well MTP	+ 6 Vials
		2 \times 48 PCR strips	+ 6 Vials
		4 \times 24 PCR strips	+ 6 Vials
Autosampler cooling	20 $^{\circ}$ C below ambient, typical setting: 7 $^{\circ}$ C (45 $^{\circ}$ F)		

Technical Specifications

Technical specifications are subdivided as follows:

- General specifications, [Table 16](#)
- Physical specifications, [Table 17](#)
- Electrical specifications, [Table 18](#)
- Communication specifications, [Table 19](#)
- Parts list, [Table 20](#)

Table 16. General

Item	Specification
Sound pressure level	< 70 dBA
Operating temperature	5 to 30 $^{\circ}$ C
Storage temperature	-25 to +60 $^{\circ}$ C
Humidity	20 to 80% RH, non-condensing
Sample viscosity	0.1 to 5 cP
Solvent A	Water with 0.1% formic acid
Solvent B	Acetonitrile with 0.1% formic acid
Safety	According to IEC 61010

Table 17. Physical

Item	Specification
Dimensions (W×D×H)	<i>w</i> 35 cm (14 in.), <i>d</i> 38 cm (15 in.), <i>h</i> 45 cm (18 in.)
Weight	32 kg (71 lb)
Weight in the shipping container	42 kg (93 lb)

Table 18. Electrical

Item	Specification
Power requirements	120 Vac, 50/60 Hz, 250 W 230 Vac, 50/60 Hz, 250 W (For UPS dimensioning, assume 250 W)
Fuses	For 120 Vac: one T 5 AL 250 V fuse (5 × 20 mm, IEC 60127) For 230 Vac: one T 2.5 AL 250 V fuse All fuses are UL Listed and CSA-certified.

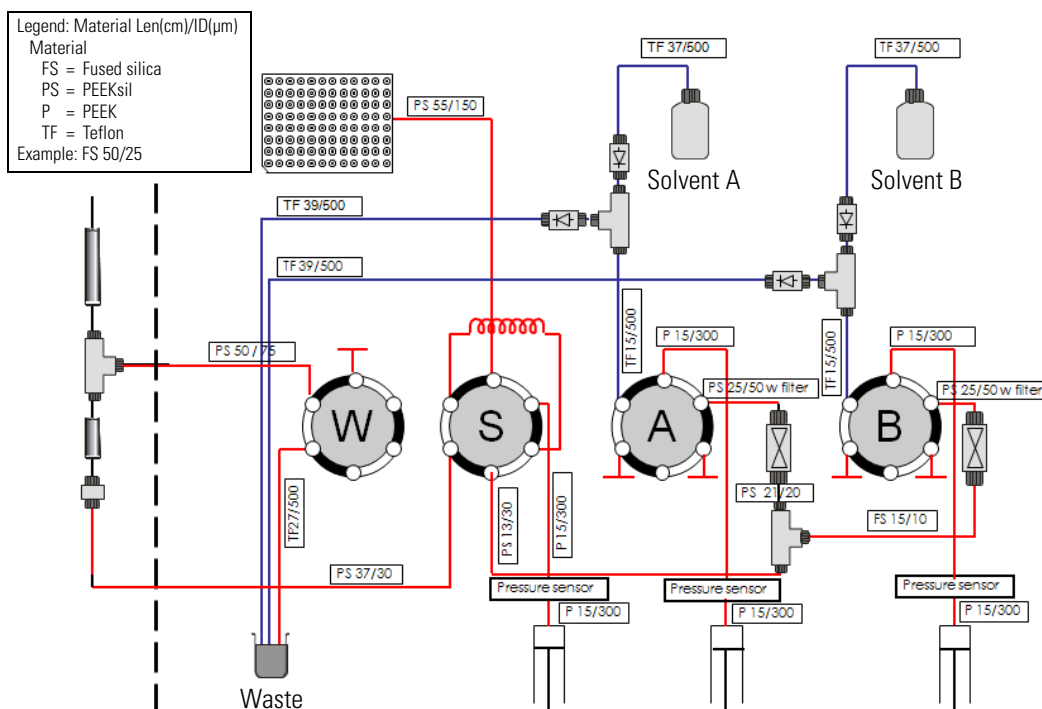
Table 19. Communication

Item	Specification
Contact closure	3 outputs, 3 inputs, and 6 ground pins
IN circuit	TTL Level
OUT circuit	PhotoMOS™ relay protected against high voltages. Continuous switch current from 0.8 to 5 V is 35 mA.
LAN	10/100 Mb/s BaseT Ethernet
USB	2 × USB 1.1 for keyboard and mouse
RS-232	Reserved for HPLC add-ons or high-level MS control by using serial communication
P-Bus	Reserved for HPLC add-ons, using the internal EASY-nLC instrument protocol bus: 8-wire control and limited power at 9/24V
PS/2	Input for connection of keyboard or mouse
Monitor	Output for connection of external display

Table 20. Parts

Item	Specification
Pumps	<ul style="list-style-type: none"> • 140 μL volume (enough for a >10 h, 0–100% B, 300 nl/min gradient) • 1 nL/min to 533 μL/min flow range • External pressure sensor
Valves	<ul style="list-style-type: none"> • VICI™/Valco™ rotor/stator • 6 ports • 3 positions (1-6, 1-2, or CENTERED)
Autosampler	<ul style="list-style-type: none"> • Peltier-cooled. Capacity is maximum 20 °C below ambient temperature, measured at 60% RH. Expect less if you remove the side panels, use an adapter plate in the autosampler other than the standard plate, or both. • Plate holder ejects through spring-mounted autosampler door. • 4 containers for waste or wash liquids

Flow Line Diagram



Miscellaneous

This appendix contains information you might refer to periodically, including how to replace the main power fuse, how to install an external device such as Advion RePlay™, a consumable parts list, and transport and disposal instructions.

Contents

- [Cleaning Materials](#)
- [EASY-nLC II Consumable Parts List](#)
- [Replacing the Main Power Fuse](#)
- [Installing and Using the RePlay External Device](#)
- [Declaration of Contamination](#)
- [Transport Instructions](#)

Cleaning Materials

When cleaning the outside of the EASY-nLC II instrument, use a mild detergent and a clean cloth.

EASY-nLC II Consumable Parts List

These EASY-nLC II parts are consumable parts and not covered by the normal first year Limited Warranty offered by Thermo Fisher Scientific, or any other service contract agreement containing extended warranty coverage.

- All flow lines including Tee-pieces and fittings (nuts, ferrules, sleeves, and valve blank stops)
- Check valves
- Solvent filters, inline filters, and all associated filter holders
- Pump piston seals
- Valve rotors and stators

- All bottles and lids
- All sample vials, microtiter plates, and associated lids and mats
- Columns
- Fuses

For consumable and spare parts descriptions and ordering information, go to www.proxeon.com/productrange/nano_lc/spares_accessories.

Replacing the Main Power Fuse

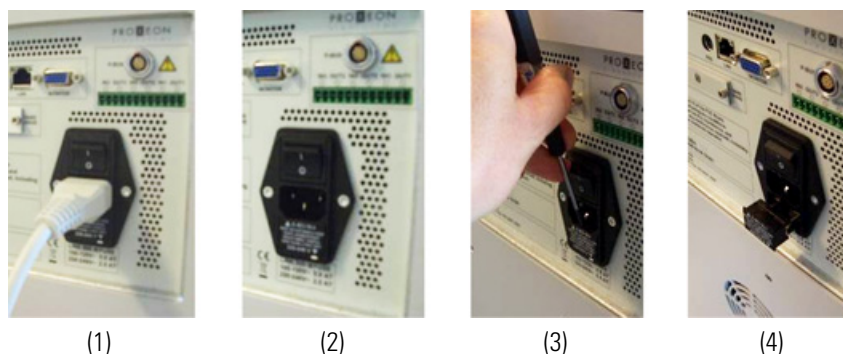
The first four numbered steps of this procedure correspond to the series of pictures in [Figure 65](#).

Note Before proceeding, make sure to turn off the system and remove the power cable.

❖ To replace the main power fuse

1. Place the EASY-nLC II instrument so that you can access the connection panel on the back side.
2. Pull out the power plug.
3. Place a screwdriver in the top opening of the fuse holder.
4. Pull the fuse holder out.

Figure 65. Replacing the main power fuse



5. Replace with the right type of fuse (see [Table 18](#) on [page 121](#)).

You can insert the fuse holder two ways:

- With the 110–120V showing at the bottom
- or–
- With the 220–240V showing at the bottom ([Figure 66](#))

Figure 66. Positioning of the 220–240V fuse



Installing and Using the RePlay External Device

Currently, the EASY-nLC II system supports the Advion RePlay™ device. Thermo Fisher Scientific may choose to support additional external devices for use with the EASY-nLC II instrument in future releases.

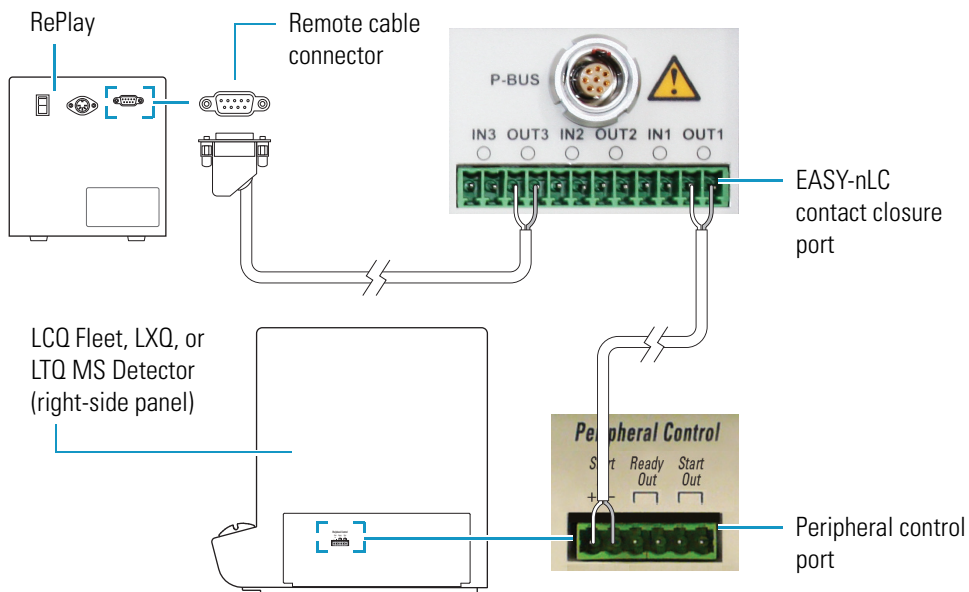
The RePlay device was specifically developed for nano LC/MS analysis. It splits the sample from one injection, so a second analysis is provided with no additional sample required. RePlay is composed of a primary column, a 6-port valve, a flow sensor, tubing, a control panel, and a secondary column/emitter. It performs a second analysis of one sample without a reinjection, which allows for twice the analytical possibilities from every nanoLC injection. For detailed information, visit the Advion BioSciences Web site:

www.advion.com/biosystems/replay/index.php

Disclaimer: Thermo Fisher Scientific does not offer support for—or is in any way responsible for—the functionality of the RePlay device itself and the associated tubing and columns that it requires.

You must connect the RePlay device to the EASY-nLC II contact closure port by using the OUT3 pins on the back panel (see [Figure 67](#)).

Figure 67. RePlay device connection to the EASY-nLC contact closure port

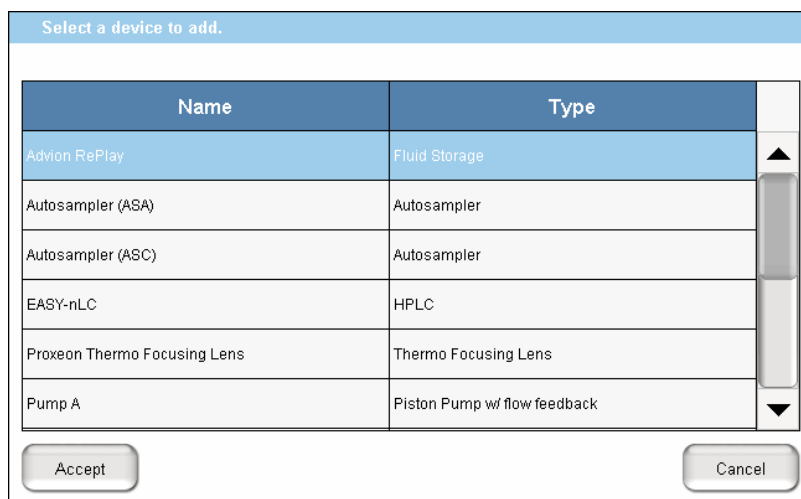


❖ **To connect RePlay to the EASY-nLC II system**

1. Turn off the instrument.
2. Connect an EASY-nLC II analytical column to RePlay valve position 5.
3. Connect Valve (4) to the RePlay focusing column and MS emitter.
4. Turn on the instrument.

❖ **To configure the RePlay device in the EASY-nLC II application**

1. Choose **Maintenance > Devices** and check that the RePlay device is not already present in the Devices list. If not, press **Add Device**.
2. In the Select a Device to Add dialog box, locate the Advion RePlay device.



3. Select the device and press **Accept**.
4. Verify that the RePlay device was added to the system by locating it in the Devices list. Selecting the device shows information about the device driver (see [Figure 68](#)).

Figure 68. RePlay About tab under the Devices list

The screenshot displays the Thermo Scientific software interface. At the top, there are navigation tabs: Home, Batch Setup, Method Editor, Maintenance (selected), and Configuration. On the left side, there is a sidebar menu with options: Scripts, Log Book, Support, and Devices (selected). Below the sidebar, there is a 'Job' section showing '00:00:09' and a date/time stamp '25 aug 2010 09:11' with the user 'admin' and an 'Exit...' button.

The main area is titled 'Devices' and contains a table with the following data:

Name	Type
Piston Pump [Pump S]	Piston Pump
RePlay	Fluid Storage
Rotary Valve [Valve A]	Rotary Valve

To the right of the table are three buttons: 'Add Device', 'Remove Device', and 'Assign Slave'. Below the table is an 'About' tab for the selected 'RePlay' device, which displays the following information:

RePlay

- Manufacturer:** Advion BioSciences, Inc.
- Product:** RePlay
- Type:** Fluid Storage
- Driver info:** 0.1.0
- Firmware info:** <N/A>
- Serial Number:** <N/A>
- Usages:** Capture/Playback Eluent
- Slaves:** <N/A>

B Miscellaneous

Installing and Using the RePlay External Device

❖ To set up a method with RePlay

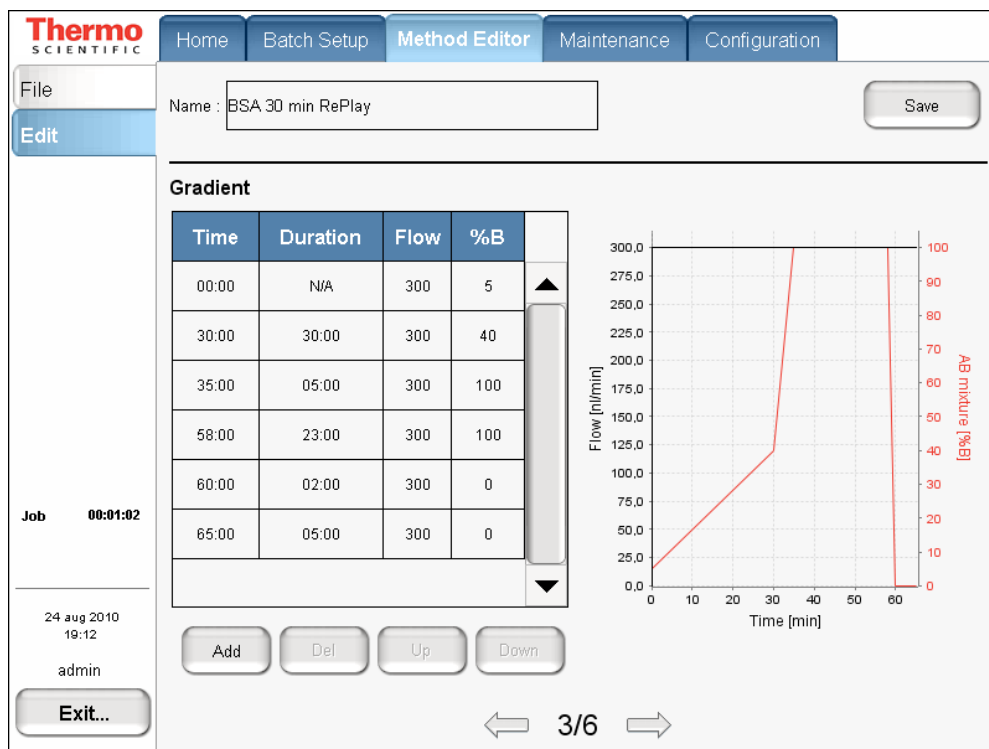
1. To use the RePlay device, enable the method extension on page 1 of the Method Editor (you can also open an existing method and enable its extension). An extra RePlay Method Editor page should now be available.

The screenshot displays the Thermo Scientific Method Editor interface. At the top, there are navigation tabs: Home, Batch Setup, Method Editor (selected), Maintenance, and Configuration. The main area is divided into several sections:

- File:** A text input field contains "Name : BSA 30 min". To its right are buttons for "Open", "Close", and "Save".
- Edit:** A blue button on the left side of the main area.
- Description:** A text area containing "Standard 30 min BSA separation w/RePlay".
- Method extensions:** A section with a checked checkbox for "RePlay" and a radio button selected for "Capture/Playback Eluent".
- Job:** A section showing "Job 00:01:02".
- Footer:** A section showing the date and time "24 aug 2010 18:34", the user "admin", and an "Exit..." button.

At the bottom right of the interface, there are navigation arrows and the page indicator "1/6".

2. On page 3 of the Method Editor, properly set up the gradient. For more information, see “To build a gradient” on page 57.



RePlay uses a nanoflow splitter and a valve to split the EASY-nLC II eluent to a direct online analysis and to a Capture Cartridge in the RePlay device. The RePlay valve switches to direct the captured chromatogram to the MS for a second analysis of the same injection. To acquire the second analysis, the gradient has to be prolonged.

3. On page 6, enter the playback delay relative to the MS START signal (which can be the start of either sample loading onto the columns or the start of the gradient, and set from the Configuration > Connections page). This value, as shown in [Figure 69](#), indicates how long into the gradient to switch back to playback mode. At the end of the gradient, the contact closure control reverts to capture mode.

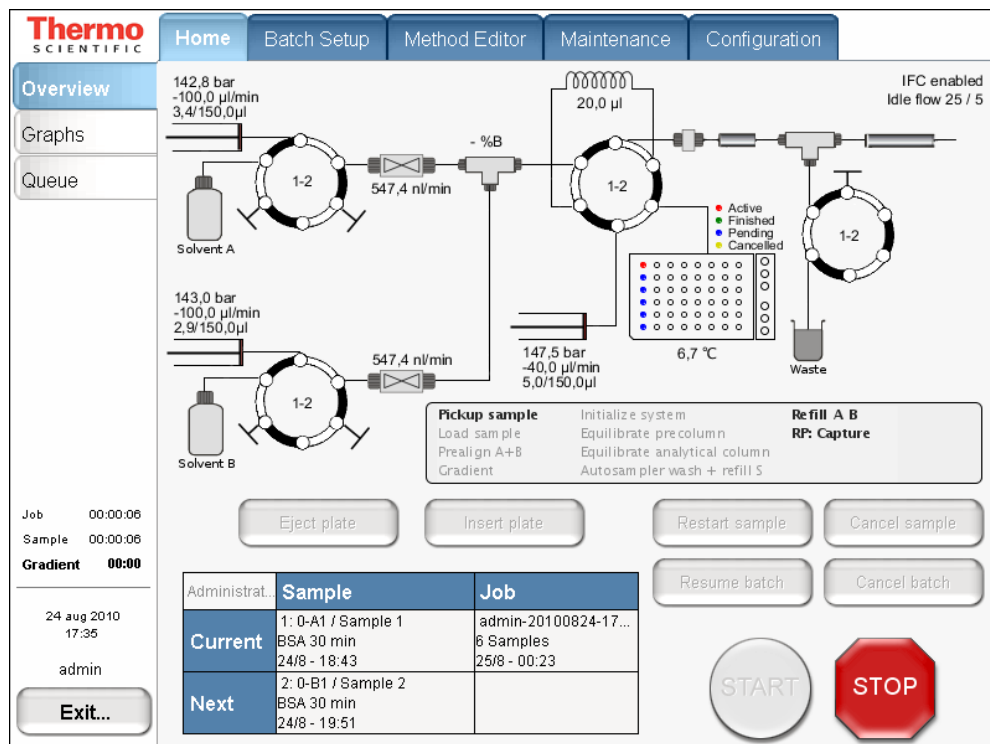
Figure 69. RePlay Playback Delay setting

The screenshot displays the Thermo Scientific software interface. At the top, there are navigation tabs: Home, Batch Setup, Method Editor (selected), Maintenance, and Configuration. Below the tabs, there is a 'File' menu and an 'Edit' button. The main area is titled 'RePlay (Capture/Playback Eluent)'. Under the 'Timing' section, the 'Playback delay' is set to '35:00 min:sec'. A 'Save' button is located in the top right corner. On the left side, there is a 'Job' status showing '00:01:02', a date and time '24 aug 2010 19:13', and the user 'admin'. At the bottom left is an 'Exit...' button, and at the bottom right are navigation arrows and the page number '6/6'.

❖ **To monitor the run of a RePlay method**

1. When running a sample using a method with the RePlay extension, open the **Overview** page under Home.
2. View the RePlay (“RP”) mode display in the center status box (see [Figure 70](#)).

Figure 70. RePlay method status display



Overhead of Using the EASY-nLC II System to Control RePlay

For a typical RePlay application—injecting and loading 5 µl of sample with a gradient of 60 minutes + 30 minutes RePlay—the full timing is as follows:

- 2 minutes – Refilling pumps at 100 µl/min
- 2 minutes – Injecting 5 µl at 20 µl/min and waiting for pressures to settle
- 12 minutes – Loading the sample onto the analytical column with 15 µl at 1.25 µl/min
- 60 minutes – Gradient + autosampler wash
- 30 minutes – RePlay while also reequilibrating the analytical column

In total, you can achieve an MS utilization rate of close to 85 percent (90/106), despite the fact that the EASY-nLC II system cannot overlap (parts of) two runs in parallel.

B Miscellaneous

Installing and Using the RePlay External Device

Column switching goes one step further in efficiency, but at the cost of extra pumps, columns, and tubing, which increases the price of the solution and introduces extra dead-volume and delays into the transfer lines.

Working with nano-flow systems, you must minimize dead-volume wherever possible and limit connections and switching as much as possible.

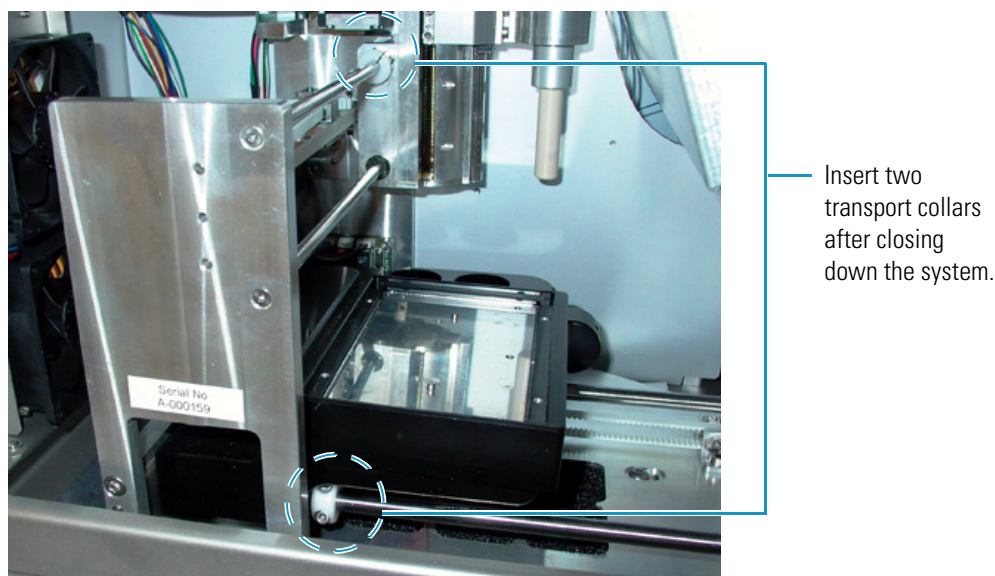
Transport Instructions

❖ To prepare for transport

1. Open the tray door on the Home > Overview page. Press **Eject Plate**.
2. Remove the plates/tubes.
3. Remove the waste container and other bottles (see “Autosampler Bottles” on page 44).
4. Close the tray door on the Home > Overview page. Press **Insert Plate**.
5. Close down the system properly (see “Closing Down the EASY-nLC II System” on page 33).
6. Remove the cables on the back side.

Tip For the original ASA autosampler only: Secure the autosampler by using the transport collars (see Figure 71) before shipping the instrument. (This is not necessary on the latest ASC model.)

Figure 71. (ASA autosampler only) Transport collar location



7. Fill out the Declaration of Contamination of Equipment in “Declaration of Contamination” on page 133.
8. Place the EASY-nLC II instrument in the original shipping container.

Index

A

accounts, adding new 40
adapter plate, replacing
 ASA autosampler 72
 ASC autosampler 74
admin, changing password 39
advanced flow control system 5
AFC *See* pump flow control
analytical Col Equilibration script 109
analytical column
 equilibrating 52
 installing 50
autosampler
 bottles, preparing 44
 calibration 71
 checking pickup 104
 setting to pick up sample 55
 Torque script 117
 troubleshooting 90
 wash 59

B

Back Pressure script 116
Batch Setup tab 23
batch, troubleshooting 88
building a gradient 57

C

calibrating
 autosampler 71
 plates 78
 wash/waste/vial positions 83
change administrator password 39
checking
 sample pickup 84
 solvent levels 103
 valves 104
column wash 60
columns, installing 50

compliance, electromagnetic *xiv*
components
 system 2
 troubleshooting 93
Configuration tab 29
configuring
 devices 40
 EASY-nLC 32
 RePlay 126
connecting
 add-ons 14
 mouse and keyboard 14
 the mass spectrometer 35
 to the network 36
connection instructions 13
consumable parts list 123
contact closure, troubleshooting 92
control system
 advanced 5
 Intelligent 6
cooler temperature, checking 104
creating
 a batch with the Microtiter plate 60
 a method 54
 new user accounts 40
 plate formats 75

D

daily maintenance 103
data types graphs 65
Declaration of Conformity *xiv*
Declaration of Contamination 133
devices, configuring 40

E

EASY-nLC II description 2
equilibrating the analytical column 52
external connections 13

F

flow control system 5
flow line diagram 122
flow parts and lines 15
flow sensor calibration 104
Flow Sensors script 117
Flush Air script 107
fuse, replacing 124

G

gradient, building 57
graphical data types 65

H

hardware overview 7
Home tab 22

I

IFC *See* pump flow control
installing columns 50
instrument
 back pressure, checking 104
 operation 4
Intelligent flow control system 6
interface layout 20
Isocratic Flow script 110

K

keyboard connection 14

L

Leaks script 113
leaks, checking for 104
lifting instructions 12
low pressure solvent filters 104

M

maintenance
 daily 103
 quarterly 104
 scripts 46, 105
 yearly 104
Maintenance tab 27
mass spectrometer connection 35
menu structure 21
Method Editor tab 25
method, creating 54

Microtiter plate 60
monitoring the run 63
 See also RePlay monitor method
mouse connection 14
MS Connection script 111

N

network access, troubleshooting 92
network connection 36

P

performance specifications 119
picking up a sample 55
plate formats 75
plates, calibrating 78
power requirements 11
pre-column equilibration 58
Precolumn Equilibration script 108
pre-column preparation 51
preparing
 autosampler bottles 44
 pre-column 51
 solvent bottles and waste containers 43
 waste container 45
 well plate 54
product description 2
pump flow control
 advanced 5
 description 5
 Intelligent 6
Purge Solvent script 106
purging and flushing pumps 47

Q

quarterly maintenance 104

R

replacing main power fuse 124
RePlay
 configuring 126
 connecting to EASY-nLC 125
 monitor method 131
 overhead of using 131
 setup method 128
RePlay (external device), installing and using 125
RePlay method 131
reset pressure sensor script 118
running batch, editing 67

S

- safety standards (EMC) [xiv](#)
- sample
 - acquisition [63](#)
 - loading, setup [55](#)
 - pick-up check [84](#)
 - Pickup script [112](#)
 - run, troubleshooting [70](#)
- saving system files on removable storage [99](#)
- scripts
 - Analytical Col Equilibration [109](#)
 - Autosampler Torque [117](#)
 - Back Pressure [116](#)
 - Flow Sensors [117](#)
 - Flush Air [107](#)
 - Isocratic Flow [110](#)
 - Leaks [113](#)
 - Maintenance [105](#)
 - MS Connection [111](#)
 - Precolumn Equilibration [108](#)
 - Purge Solvent [106](#)
 - Reset Pressure Sensor [118](#)
 - Sample Pickup [112](#)
 - Valve Check [114](#)
 - Valve Tune [117](#)
- setup
 - mass spectrometer connection [35](#)
 - method with RePlay [128](#)
 - pre-column equilibration [58](#)
- solvent bottles, preparing [43](#)
- solvent levels, checking [103](#)
- specifications
 - performance [119](#)
 - technical [120](#)
- starting EASY-nLC [19](#)
- stopping sample acquisition [66](#)
- system configuration [3](#)
- system files, saving to removable storage [99](#)
- system overview, monitoring [64](#)
- system startup, troubleshooting [93](#)

T

- technical specifications [120](#)
- temperature and humidity [12](#)
- touch screen
 - interaction principles [21](#)
 - using [18](#)
- transport instructions [134](#)
- troubleshooting
 - autosampler [90](#)
 - contact closure [92](#)

- network access [92](#)
- running a batch [88](#)
- sample run [70](#)
- system components [93](#)
- system startup [93](#)
- turning off EASY-nLC [33](#)

U

- user interface layout [20](#)
- user permissions [38](#)

V

- Valve Check script [114](#)
- Valve Tune script [117](#)
- verifying mass spectrometer connection [63](#)

W

- warranty [123](#)
- wash column [60](#)
- waste container, preparing [43, 45](#)
- well plate, preparing [54](#)

