

Installation Manual for AAA Direct

023741 Revision 05 • July 2018



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

Installation Manual

for

AAA Direct

© 2018 Thermo Fisher Scientific Inc. All rights reserved.

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

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

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

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.

Revision History:

Revision 05, July 2018, Content reformatted to meet current branding and ISO standards. Minor corrections made to specification data.

Safety and Special Notices

Make sure you follow the precautionary statements presented in this guide. The safety and other special notices appear in boxes.

Safety and special notices include the following:



Indicates a potentially hazardous situation which, if not avoided, could result in death or serious injury.



Indicates a potentially hazardous situation which, if not avoided, could result in damage to equipment.



Indicates a potentially hazardous situation which, if not avoided, may result in minor or moderate injury. Also used to identify a situation or practice that may seriously damage the instrument, but will not cause injury.



Indicates information of general interest.

IMPORTANT

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

Tip

Highlights helpful information that can make a task easier.

Contents

1.	Intro	duction	6			
2.	The	Most Important Rules	7			
2	2.1	Important Preparations: (Always do the following)				
2.2		Important Reminders: (Never do the following)	7			
3.	Initia	al Check List	8			
4.	Insta	llation	9			
4	.1	System Configuration and Start-up	9			
	4.1.1	System Configuration (Time Requirement: 120 min)	9			
	4.1.2	Creating a Sequence and an Instrument Method	10			
	4.1.3	System Rinse (Time Requirement: 75–90 min)	10			
4	.2	Verification of System Cleanliness	11			
	4.2.1	Eluent Preparation (Time Requirement: 45 min)	11			
	4.2.2	System Background Check (Time Requirement: 30 min)	11			
4	.3	Verification of System Response	11			
	4.3.1	Adjusting Eluent Composition (Time Requirement: 10 min)	11			
	4.3.2	Column Installation (Time Requirement: 15 min)	12			
	4.3.3	Histidine Injection (Time Requirement: 30 min)	12			
4	.4	Verification of System Functionality	14			
	4.4.1	Injection of NIST SRM 2389 Standard (Time Requirement: 100 min)	14			
5.	Troul	bleshooting	16			
5	5.1	Verification of System Cleanliness	16			
5	5.2	Verification of System Background	16			
5	5.3	Verification of System Functionality	17			

1. Introduction

Amino acid analyses require attention to detail, regardless of the analytical approach. The Dionex AAA-Direct system simplifies that task because it eliminates the need for additional high purity reagents that are required by other techniques.

This procedure is written specifically for Dionex service representatives working closely with the nearest Dionex Applications Lab to ensure a successful installation experience. Please read the entire document carefully before arranging an installation.

2. The Most Important Rules

2.1 Important Preparations: (Always do the following)



Wear appropriate gloves (non-powder) when handling eluent bottles, samples, or electrode cell parts. Don't touch these with your bare hands. Gloves will protect you from the chemicals and protect the parts from fingerprint damage.

- Use 50% NaOH solution rather than NaOH pellets to make eluent.
- Use dedicated glassware and disposable glass or plastic ware for volume adjustments.
- Keep your NaOH eluent blanketed by inert gas. Prepare new NaOH eluent if left unblanketed for more than 30minutes.
- Prime at least 40 mL of new eluent through the lines when changing eluent. This will ensure that your fresh eluent is primed through the lines up to the pump heads.
- Use pre-slit septa with the injection vials.
- Use 25 µL loop size; larger loops will cause loss of resolution.
- Install and use the piston wash option.

2.2 Important Reminders: (Never do the following)

- Do not continue to the next step of the procedure if the previous has failed.
- Do not start an installation with any of the checklist items missing (see Section 3 for Check List).

Never use bottled HPLC water. Do not store deionized water (DI water), always use freshly drawn DI water for any preparation of eluents.

Never use 'communal' filtration units or filters made of unknown or unsuitable (cellulose derivatives, polysulfone) materials.

• Never use eluent inlet filters; cover the ends of the eluent lines with parafilm when changing bottles.

Never use MeOH or other organic solvent as rinse fluid in the autosampler. Use only 20 ppm sodium azide, or water if replaced weekly.

• Do not run above 50 °C or 3,500 psi.

3. Initial Check List

These items *MUST* be available at the customer site. Check with the customer when you call to schedule the installation.

- □ Laboratory water unit delivering 18.2 megohm-cm water at the installation site.
- □ Vacuum pump available for use with the vacuum filtration units.
- □ Three sterile-packed Nylon Nalgene Filtration Units, Funnel Size 1.0 L (VWR Item # 28198-514).
- \Box Inert gas cylinder (helium or nitrogen) with a regulator valve (about 0–200 psi at the low-pressure side) and the appropriate size adaptors plus tubing.
- □ Dionex AAA-Direct Installation Kit (Item # 059539) containing: This Standard Operating Procedure, Histidine quality solution, two bottles of sodium acetate, one bottle of 50% sodium hydroxide, and restriction tubing.
- □ Product Manual for the AAA-Direct (Document. No. 031481) available on https://www.thermofisher.com/search/results?query=aaa%20installation%20manual&focu sarea=Search%20All.
- □ NIST Amino Acid standards (SRM 2389, 2.5 mM solution).
- \Box One spare AAA certified Au electrode (Item # 063722).
- \Box One spare pH-Ag/AgCl reference electrode Item # 044198).
- Sterile-packed, 10 mL and 25 mL disposable pipets and suitable pipetting bulbs or pumps.
- □ Sodium azide solid, reagent grade for preparation of diluent solution.
- \Box Powder-free, disposable gloves (at least 1 box).
- Disposable, plastic (PE) large-size (at least 20 mL) syringe for priming the pump.
- □ Nitric acid, 2 N, concentrated nitric acid, diluted 1:1 (v/v) for system cleanup.

4. Installation

Make sure that each of the following sections are completed successfully before moving onto the next section. If you are having problems, check the troubleshooting guide at the end of this procedure. If you are still having problems, call Dionex.



Refer to the AAA-Direct manual for assistance and precautions.



The instructions in this manual are for use with non-disposable gold working electrodes only (Item # 063722). Please refer to the AAA-Direct Product Manual (Document No. 031481) when working with disposable electrodes (Item # 060082 - 6 electrode and Item # 060140 - 24 electrode).

4.1 System Configuration and Start-up

4.1.1 System Configuration (Time Requirement: 120 min)



Never polish a disposable gold electrode.

- a) Configure the system as follows:
 - 1. The AS-AP or equivalent autosampler on the top of the DC
 - 2. The pump on the right (on the top of EG module if available).
- b) Ensure that nitrogen or helium is delivered to the eluent organizer with about 5–6 psi at each bottle.
- c) Check that the AS-AP or equivalent is plumbed with red tubing (0.005 i.d).
- d) Check that the red tubing is minimized to reduce dead volume.



Do not connect the column as that will be completed in a future step.

- e) Make all fluid and electrical connections.
- f) Install the yellow tubing from the Installation Kit between the injector and detector cell inlet.
- g) Assemble the electrochemical cell with the gold (Au) AAA-Direct-Certified working electrode (*Item # 063722*).
- h) Verify that the modules are communicating.

4.1.2 Creating a Sequence and an Instrument Method

Create a sequence "HisNIST" and two Instrument Methods by following the example in Appendix B. The waveform can be programed into the Instrument Method by selecting "pH" as "Reference Electrode" and "Gold, pH-Ag-AgCl, RE, AAA" as "Waveform" in the "EDet1 Options for Options (ED:EDet1)" during the creation of Instrument Method in the Chromeleon (CM) software. Please refer to the manual of CM for more detail about it (Quick Start Guide Chromeleon 7.2, Item #: 7, rev 1.0, 2013). The waveform is also shown in Table 1 (page 26) and the gradient profile is from table 4 (page 49) in Product Manual for AAA-Direct (031481-14, 2018).

Figure 1 HisNIST Sequence as it should appear in the Data Directory following Restore

Sequen	ce: HisNIST	HisNIST							Page 1 of 1	
Operato	or: jcheng	jcheng Print							ed 7/2/01 5:15:43 PM	
Title: Datasou Locatio Timeba #Sampl	eds seque urce: PJANDI n: sequence sse: PJANDI les: 4	ence K_SYS2 \Jul2001\Car K_SYS2_1	eKit		Created: Last Upd	ate:	6/28/01 5:48:10 PM by jcheng 7/2/01 1:37:53 PM by jcheng			
No.	Name	Type	Pos	Inj.	Program	Method	l Status	Inj. Date/Time	Comment	
1	WaterBlankIsocratic	Unknown	1	25.0	36A24B40C_10min	His	Finished	6/28/01 6:52:25 PM	Response test	
2	8uMHis	Standard	2	25.0	36A24B40C_10min	His	Finished	6/28/01 9:58:44 PM	Response test	
3	WaterBlankGradient	Unknown	1	25.0	GP50Comp_010PN6	NIST	Finished	6/29/01 3:48:29 PM	System test	
4	8uMNISTsrm2389	Standard	3	25.0	GP50Comp_010PN6	NIST	Finished	6/29/01 5:04:36 PM	System test	

4.1.3 System Rinse (Time Requirement: 75–90 min)

- a) Prepare a solution of 2 M NaOH to rinse each bottle:
 - 1. Dilute 104 mL of 50% sodium hydroxide to 1 L of Deionized Water (DI water)
- b) Place the 2 M NaOH in a pre-rinsed bottle.
- c) Place all 4 eluent lines in the pre-rinsed bottle.
- d) Prime at least 40 mL of sodium hydroxide from each line.
- e) Close the solvent "draw-off" valve.
- f) Leave the pump proportioning at 25/25/25 for 15 minutes.
- g) Ensure all surfaces come into contact with the sodium hydroxide (rotate the injection valve).
- h) Repeat the process with DI Water (18.2 megohm-cm water).
- i) If everything completed successfully then continue to the next section.

4.2 Verification of System Cleanliness

4.2.1 Eluent Preparation (Time Requirement: 45 min)

- a) Prepare eluents as described in the AAA-Direct Product Manual (Document No. 031481) and transfer them into their respective eluent reservoirs making sure to follow all of the precautions.
- b) Set the eluent component to 100% for each eluent
- c) Prime at least 40 mL of eluent from each reservoir after filling the eluent bottles.

4.2.2 System Background Check (Time Requirement: 30 min)

- a) Verify system background as described in the Section 6.2.1. in the AAA-Direct Manual (031481-14, 2018).
- b) Ensure the following:
 - 1. The detector is set to pH mode (not Ag mode)
 - 2. The cell is not on
 - 3. The pump is pumping at 0.25 mL/min
 - 76% DI water (Channel A)
 - 24% 0.25 M NaOH (Channel B)
- c) Confirm the pH reading in the ED panel of the Instruments is between 12.1 and 13.0.
- d) Turn on the cell and begin monitoring the background signal from the control panel for at least 30 min.
- e) Confirm the baseline is < 80 nC. If the background exceeds 80 nC or the pH is out of range, see the Troubleshooting Section.

4.3 Verification of System Response

4.3.1 Adjusting Eluent Composition (Time Requirement: 10 min)

- a) Change eluent composition to (flow rate: 0.25 mL/min):
 - 1. 36% DI water (Channel A)
 - 2. 24% 0.25 M NaOH (Channel B)
 - 3. 40% 1.0 M NaOAc (Channel C)
- b) Wait 10 minutes until the background is stable and <130 nC.
 - If it is drifting down, wait as long as it takes to stabilize below 130 nC.
 - If the background exceeds 130 nC, see the troubleshooting section.
 - If the background is stable and <130 nC, proceed to the next section.

4.3.2 Column Installation (Time Requirement: 15 min)

- a) Stop the flow.
- b) Turn off the cell voltage.
- c) Remove the yellow restrictor tubing.
- d) Install the AminoPac PA10 guard and analytical columns, but **DO NOT** connect the column outlet to the cell inlet yet.
- e) Turn the pump back on.
- Pump the following through the column and into a waste container for 10 min at 0.25 mL/min.
 - 1. 36% DI water (Channel A)
 - 2. 24% 0.25 M NaOH (Channel B)
 - 3. 40% 1.0 M NaOAc (Channel C)
- g) Connect the column tubing to the cell.
- h) Verify the background is still <130 nC.
 - If it is not, see the troubleshooting section
 - If the background is stable and <130 nC, proceed to the next section

4.3.3 Histidine Injection (Time Requirement: 30 min)

- a) Make an 8 μ M solution of histidine by adding 1 mL of water to the dry residue in the micro vial shipped with installation kit.
- b) Place a vial with DI water in position RA1 of the autosampler.
- c) Place the histidine quality solution in positionRA2.
- d) Run lines 1 and 2 as shown in the Installation sequence created as a copy of the HisNIST sequence as described in the section 6.1.1. in the AAA-Direct Manual (031481-14, 2018).
- e) Confirm the peak height for histidine is >200 nC (see Figure 2).
- f) Ensure the % RSD for "His" peak height is <5%.
 - If this is not the case, see the troubleshooting section



Figure 2 Testing the Detection Response

4.4 Verification of System Functionality

4.4.1 Injection of NIST SRM 2389 Standard (Time Requirement: 100 min)

- a) Program the pump to deliver at 0.25 mL/min:
 - 76% DI water (Channel A)
 - 24% 0.25 M NaOH (Channel B)
- b) Let the system equilibrate.
- c) Set the column oven to 30 $^{\circ}$ C.
- d) Verify that the background level is < 80 nC.
- e) Prepare 1 L of 20 mg/L of sodium azide in DI water.
- f) Prepare 100.00 mL of 8 µM NIST standard by doing the following:
 - 1. Pipetting exactly 320.0 μL of NIST SRM 2389 concentrate into a clean 100 mL volumetric flask
 - 2. Filling up the flask to 100 mL with the 20 mg/L azide solution.
- g) Ensure that there is still a water blank (vial of DI water) in position RA1 of the autosampler.
- h) Place the 8 μ M NIST standard into position RA3.
- Execute lines 3 and 4 of the Installation sequence (please see Appendix B The Sequence and the Instrument Methods in the in the AAA-Direct Manual (031481-14, 2018).
- j) Confirm that the baseline rise from the start of the run to the top of the acetate gradient does not exceed 50 nC.
 - If this is not the case, see the troubleshooting section
- k) Confirm that the Arginine peak is >120 nC/235 pmol (see Figure 3).
 - If this is not the case, see the troubleshooting section
- 1) Overlay your separation with the information from line 4 of the HisNIST sequence.
- m) Confirm that the resolution between Ala and Thr is comparable.
 - If this is not the case, see the troubleshooting section



If everything is successful to this point then installation is complete.



5. Troubleshooting

5.1 Verification of System Cleanliness

- If pH is not within range:
- a) Confirm that the eluent composition is set to 76/24% (A/B).
- b) If eluent is correct, recalibrate Reference Electrode (see instructions in the ED50 Manual).
 - If the reference electrode cannot be calibrated, check the cable from the detector to the cell.
 - Alternatively, replace the reference electrode.
 - If background >80 nC and not decreasing:
- c) Repeat the 2 M sodium hydroxide rinse through all lines.
 - If the baseline remains at >80 nC after the second sodium hydroxide rinse, rinse the entire system, including all the bottles and eluent lines with 2 N nitric acid, overnight.
 - Replace the AAA-Direct gold working electrode with the spare if the baseline continues to exceed 80 nC.
 - If either the baseline or the pH remain out of range, call Thermo Dionex Support.

5.2 Verification of System Background

- If background remains steady but exceeds 130 nC:
- a) Re-rinse the acetate channel eluent lines and eluent bottle with 2 N NaOH.
- b) In addition, rinse the acetate channel eluent lines and eluent bottle 3-times with purified filtered water.
- c) Make up fresh acetate and try again.
 - If after the column is connected to the cell, the background exceeds 130 nC and is not dropping:
- d) Remove the column and replace the restrictor tubing, pumping 36:24:40 (A:B:C).
- e) If the background is < 130 nC, install a new AminoPac PA10.
- f) If the background exceeds 130 nC with the restriction tubing and is not dropping:
 - 1. Discard the acetate eluent
 - 2. Rinse line C and the rest of the system thoroughly with DI water from line A.
 - 3. Prepare new acetate and re-check the background.
 - If it is high again, you either have a contamination source in your labware or a bad lot of acetate.
 - Call Thermo Dionex for help.
 - If the histidine peak height is out of range:
- g) Check the injector and confirm that the 25 μ L injection loop is installed.
- h) Check if the autosampler is set properly and functioning correctly.
 - If the autosampler is set properly and functioning correctly:

- 1. Remove the column
- 2. Install the restriction tubing
- 3. Inject the histidine.
- i) The height of the histidine peak near the void should exceed 180 nC:
 - If the peak height is > 180 nC, install a new column.
 - If the peak height is < 180 nC, replace the AAA-Direct gold working electrode with the spare and repeat the installation section.

5.3 Verification of System Functionality

- If the acetate gradient causes a baseline rise exceeding 50 nC but the second injection shows less rise than the first:
- a) Run 6:24:70 (A:B:C) through the column for 1 hour
- b) Re-equilibrate for 30 minutes at 76:24 (A:B)
- c) Repeat the gradient test.
 - If the first and second injections show similar and excessive baseline rise, the eluent lines should be cleaned with 1 M NaOH as described in the AAA-Direct Manual.
 - If the arginine peak is <140 nC/200 pmol replace the AAA-Direct gold working electrode with the spare.
- d) Call Thermo Dionex should the arginine peak height be lower than 140 nC after installation of the spare Au electrode.
 - If you observe a resolution loss between Ala and Thr:
- e) Check the supply of inert gas to channel B (250 mM NaOH).
 - Carbonated NaOH will cause resolution loss and retention time fluctuation in the Ala to Met region of the chromatogram.
- f) If poor resolution persists, remove guard column and re-inject the same standard using the main column only.
- g) Replace the guard if the resolution is restored.
- h) Replace the analytical column if poor separation persists without the guard.
 - Another possible source could be the temperature accuracy of the Column Compartment of the DC. Call Dionex Thermo should poor separations persist between Ala and Thr after removing the guard.



Avoid pressure fluctuations. Severe pressure pulses can damage the guard column causing poor peak shapes (split peaks, peak fronting, tailing) and resolution.