

DIONEX GUARDCAP VIAL CAPS

QuickStart Guide

Setup of Dionex AS-DV

The default sample transfer line in the Dionex AS-DV autosampler should be replaced with a shorter and narrower reduced volume line to optimize the ion exchange capacity of Dionex Guardcap vial caps. In addition, several recommended settings should be used in the instrument method/program. Please see the tables below for more information.



NOTE

Detailed instructions for use can be found in the Dionex Guardcap Product Manual (Doc No. 065705), available online at <http://www.thermofisher.com/Guardcap>.

Recommended Reduced Volume Sample Transfer Lines

	Injection valve mounted in Dionex IC system	Injection valve mounted in Dionex AS-DV
Transfer line length (volume)	70 cm (35 µL) Part number 22120-60055	35 cm (17.5 µL) Part number 22120-60051

Recommended Instrument Method/Program Settings for Dionex AS-DV

	Injection valve mounted in Dionex IC system	Injection valve mounted in Dionex AS-DV
Sample delivery volume	350 µL	175 µL
Loading mode	Concentrator mode	Concentrator mode
Delivery speed	1.0 mL/min (0.5 mL/min for Dionex Guardcap HRP)	1.0 mL/min (0.5 mL/min for Dionex Guardcap HRP)

Filling the Dionex PolyVials and Installing Dionex Guardcap Vial Caps

1. Weigh 6.1-6.4 g sample/vial and install the Dionex Guardcap vial cap using the flat end of the installation tool so the cap is flush with the top of the vial. Note that this amount of sample is above the top of the vial tray. The backpressure applied by the user to install the Dionex Guardcap vial cap is higher than when installing a filter or plain cap. This is normal since Dionex Guardcap vial caps contain ion exchange resin in the pores of the filters.
2. After installation of the Dionex Guardcap vial cap, invert the vial to confirm that no air bubble is freely floating under the cap.
3. Before loading the vial into the carousel, shake off any liquid that has been forced into the cap socket.

Reordering Information

Part number	Description
302504	Dionex Guardcap H vial caps, 5 mL, Package of 125
302797	Dionex Guardcap Na vial caps, 5 mL, Package of 125
302799	Dionex Guardcap HRP vial caps, 5 mL, Package of 125
302982	Dionex Guardcap AEX vial caps, 5 mL, Package of 125
302765	Dionex Guardcap H vial caps, 5 mL, Package of 12 (includes 70 cm sample transfer line)
302796	Dionex Guardcap Na vial caps, 5 mL, Package of 12 (includes 70 cm sample transfer line)
302798	Dionex Guardcap HRP vial caps, 5 mL, Package of 12 (includes 70 cm sample transfer line)
302981	Dionex Guardcap AEX vial caps, 5 mL, Package of 12 (includes 70 cm sample transfer line)
22120-60055	70 cm sample transfer line for use with injection valves mounted in Dionex IC systems
22120-60051	35 cm sample transfer line for use with injection valves mounted in Dionex AS-DV