
THE TRAP COLUMN/SUPPRESSOR CLEAN-UP KIT INSTALLATION AND OPERATION MANUAL

The Trap Column/Suppressor Clean-up Kit (P/N 059659) is a stand-alone pneumatically driven reagent delivery system designed for delivery of suitable reagent solutions to regenerate and/or cleanup DIONEX IonPac Trap Columns and Suppressors. The use of this kit eliminates the need for using the analytical pump for delivery of concentrated acids, bases or other reagents that can cause serious contamination to the IC system.

The kit consists of a 1-liter plastic reservoir, a gas pressure regulator (25 psi max), a two-way liquid shut-off valve, and the necessary tubing and connection fittings; the kit is pre-assembled and easy to install and operate.

SECTION 1. INSTALLATION

1. Connect the gas pressure regulator (P/N 038201) to a pressurized helium source using a section of the gas line tubing (P/N 052295). Be sure to connect the gas line tubing to the inlet fitting of the regulator (the opposite fitting marked with an arrow on top of regulator).
2. Connect the exit side of the gas pressure regulator to the plastic bottle reservoir using a section of gas line tubing.
3. Install the end-of-line filter (P/N 045987) on the liquid line that extends into the bottle reservoir to trap any particulate matter.
4. Remove the union coupler. In its place, install the device to be treated and direct the liquid to a waste container through the green PEEK tubing. See Figure 1 "Regeneration of a Trap Column using the Trap Column/Suppressor Clean-up Kit" or Figure 2 "Clean-up of a Suppressor using the Trap Column/Suppressor Clean-up Kit" for how to make appropriate connections.

SECTION 2. OPERATION

2.1 REGENERATION OF IONPAC TRAP COLUMNS

1. Prepare the correct regenerant solution according to Document No. 032697 "Installation Instructions and Troubleshooting Guide for the IonPac Anion Trap Columns" or Document No. 034536 "Installation Instructions and Troubleshooting Guide for the IonPac Cation Trap Columns" or Document No. 034990 "Installation Instructions and Troubleshooting Guide for the IonPac MFC-1 Metal-Free Trap Column." Place the regenerant solution in the plastic reservoir and pressurize the bottle using the pressure regulator.
2. Direct the liquid line to a waste container and open the shut-off valve. Adjust the pressure regulator to 10-15 psi for most cases. If necessary, the pressure may be increased to a maximum of 25 psi. A safety relief valve on the regulator prevents pressure greater than 25 psi from being applied to the regenerant reservoir.
3. Deliver the recommended total volume of regenerant solution through the trap column. Refer to the corresponding product manual for the total volume recommended for the given trap column.
4. When completed, remove the regenerated trap column and install it back in the IC system. Follow the steps described in the corresponding trap column QuickStart and/or product manual to re-start the IC system.

2.2 CLEAN UP OF SUPPRESSORS

1. Refer to the suppressor clean-up section of the corresponding suppressor product manual for the correct contamination clean-up procedure.
2. Prepare and place the clean-up solution in the plastic reservoir and pressurize the bottle using the pressure regulator.
3. Direct the liquid line to a waste container and open the shut-off valve. Adjust the pressure regulator to 15-20 psi for most cases. If necessary, the pressure may be increased to a maximum of 25 psi. A safety relief valve on the regulator prevents pressure greater than 25 psi from being applied to the regenerant reservoir.
4. Deliver the recommended cleaning solution through the suppressor. Refer to the clean-up procedure in the corresponding product manual for the recommended amount of solution for the given type of suppressor.
5. When the clean-up is completed, rinse the suppressor with deionized water. Follow the steps described in the clean-up section of the corresponding product manual to start the IC system.
6. Perform the recommended QuickStart procedure for the suppressor before installation in the IC system.
7. Install the suppressor and restart the IC system.

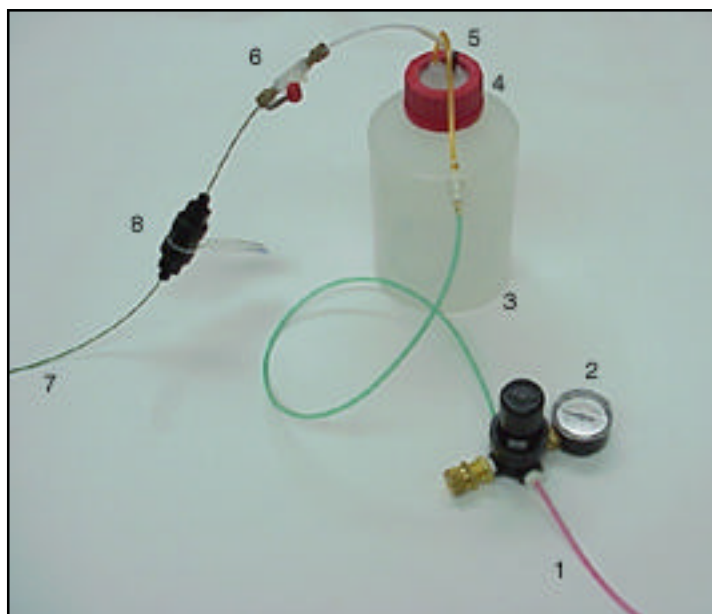


Figure 1

Regeneration of a Trap Column Using the Trap Column/Suppressor Clean-up Kit

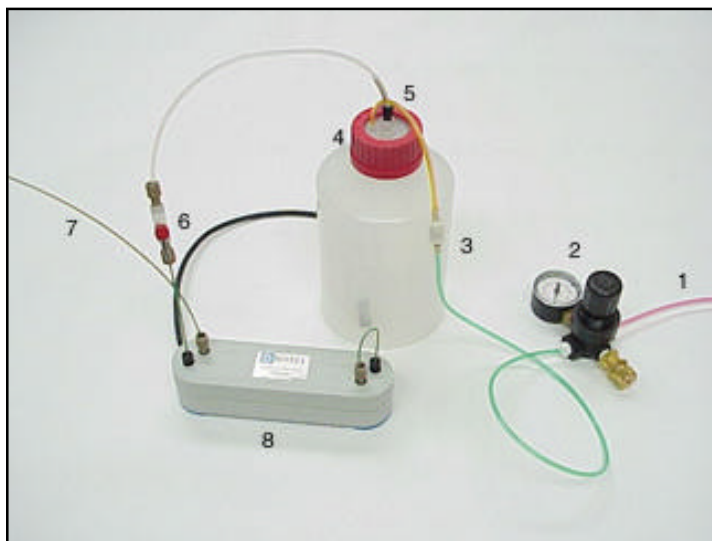


Figure 2

Clean-up of a Suppressor Using the Trap Column/Suppressor Clean-up Kit

Item	Description	P/N
1.	Air line tubing, 10 ft	052295
2.	Pressure regulator, 25 psi	038201
3.	Plastic bottle, 1liter	045902
4.	Bottle cap with bore	045897
5.	Bottle stopper	057715
6.	2-way valve	059650
7.	PEEK tubing, 0.030" I.D., 5 ft	052305
8.	Suppressor or Trap Column	
	End of line filter	045987 Not shown
	PEEK Tube cutter	045257 Not shown
	Replacement blades for PEEK Tube cutter	045258 Not shown

Now sold under the
Thermo Scientific brand

Thermo
SCIENTIFIC

Part of Thermo Fisher Scientific