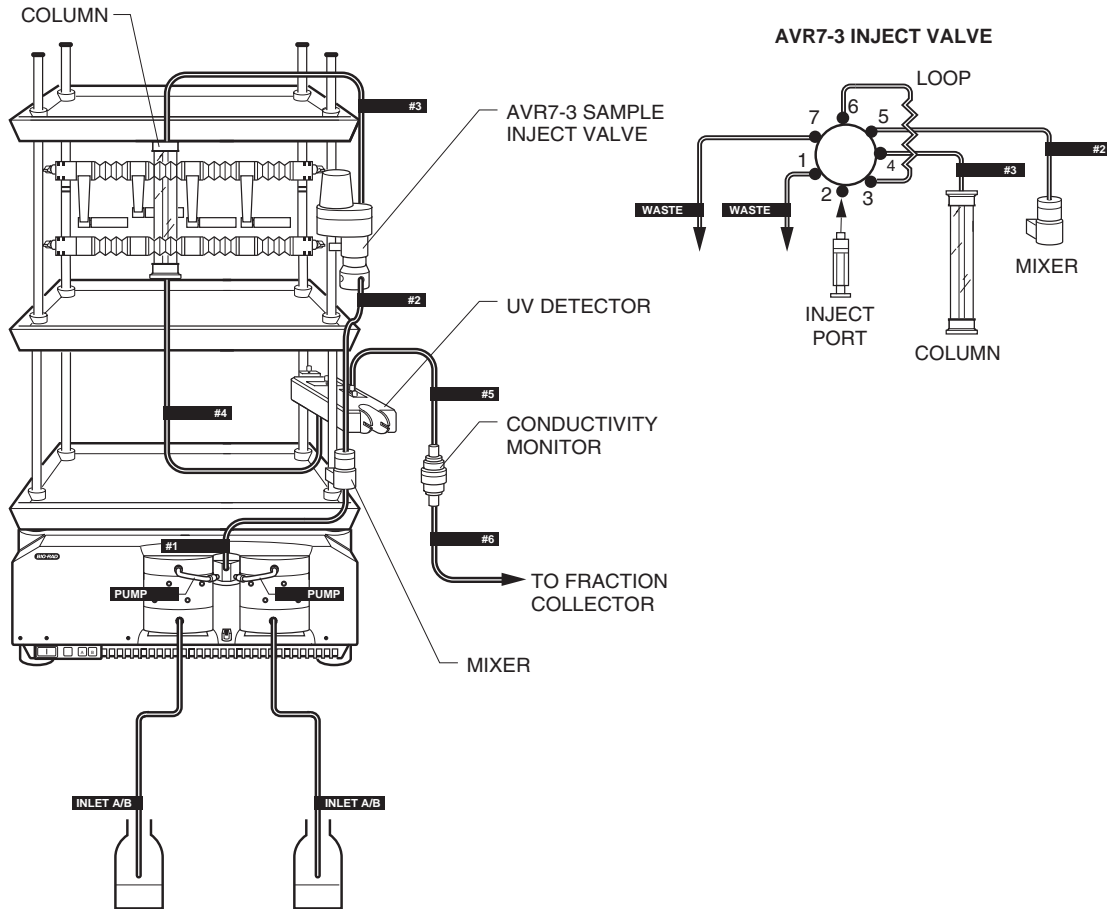


BioLogic DuoFlow F40 Tubing Kit

Catalog Number 760-0652

Dear Customer,
 The BioLogic DuoFlow F40 tubing kit is a set of pre-made tubing to simplify initial setup of your system when a large column that produces high back pressure is in use. The recommended maximum back pressure for this tubing is 3000 psi.

Follow the diagram below to complete the installation. Note that orange PEEK tubing identical to that labeled "Pump" in this kit may already be installed on the Workstation. Depending upon the flow rates and amount of back pressure produced by your system orange PEEK may still be desirable between the sample injection, column, and detector.



Tubing Label	Tubing Location	Quantity x Tubing Length	Tubing Material, OD, ID	Fittings
Inlet A/B	Buffer containers to pump inlets A and B	2 x 100 cm ± 1 cm	PTFE, clear OD=1/8" ID = 0.062"	¼ -28 black Delrin fitting on one end of the tubing only.
Waste	From ports #1 and #7 of the AVR7-3 valve to waste container	2 x 100 cm ± 1 cm	Tefzel, clear OD = 1/16" ID = 0.020"	¼ -28 gray Delrin fitting on one end of the tubing only.
Pump	From pump outlets to pressure transducer inlets (note: this tubing may already be installed on your Workstation)	2 x 10 cm ± 1 cm	PEEK (Green) OD = 1/16" ID= 0.030"	¼ -28 gray Delrin fittings on both ends of the tubing.
# 1	From pressure transducer outlet to dynamic mixer inlet	1 x 28 cm ± 10 cm	PEEK (Green) OD = 1/16" ID= 0.030"	¼ -28 gray Delrin fittings on both ends of the tubing.
# 2	From dynamic mixer outlet to port 5 of AVR7-3 valve	1 x 360 mm ± 1 cm	PEEK (Green) OD = 1/16" ID= 0.030"	¼ -28 gray Delrin fittings on both ends of the tubing.
# 3	From port #4 of the AVR7-3 valve to column inlet	1 x 24 cm ± 1 cm	PEEK (Green) OD = 1/16" ID= 0.030"	¼ -28 gray Delrin fittings on both ends of the tubing.
# 4	From column outlet to UV detector inlet	1 x 24 cm ± 1 cm	PEEK (Green) OD = 1/16" ID= 0.030"	¼ -28 gray Delrin fittings on both ends of the tubing.
# 5	From UV detector outlet to conductivity monitor inlet	1 x 15 cm ± 1 cm	PEEK (Green) OD = 1/16" ID= 0.030"	¼ -28 gray Delrin fittings on both ends of the tubing.
# 6	From conductivity monitor to the fraction collector	1 x 52 cm ± 1 cm	Tefzel, clear OD = 1/16" ID = 0.020"	¼ -28 gray Delrin fitting on one end of the tubing only.