



## BioLogic DuoFlow™ F10 Tubing Kit

Catalog #760-0650

The BioLogic DuoFlow F10 tubing kit comprises premade tubing to simplify initial setup of your system. The recommended maximum backpressure for this tubing is 5,000 psi.

Follow the diagram below to complete the installation.

**Note:** Tubing identical to that labeled **Pump** in this kit may already be installed on the workstation.

### BioLogic DuoFlow chromatography system

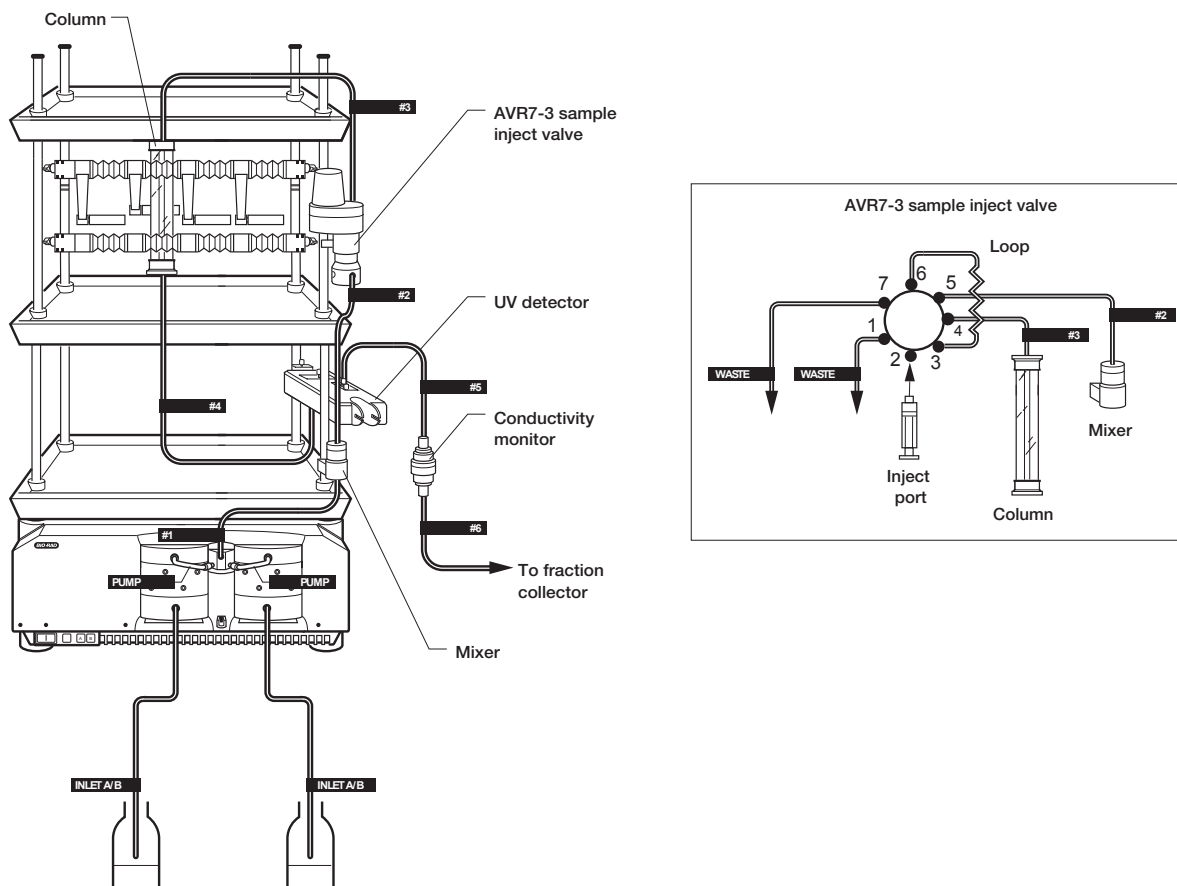


Fig. 1. BioLogic DuoFlow chromatography system tubes. Diagram shows how each labeled tube from the kit is plumbed. Inset: AVR7-3 sample inject valve, enlarged.

**Table 1. Tubing label and location.**

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"	1/4-28 black Delrin fitting on one end of the tubing only
Waste	From ports #1 and #7 of AVR7-3 valve to waste container	2 x 100 cm ± 1 cm	Tefzel, clear OD = 1/16" ID = 0.020"	1/4-28 black Delrin fitting on one end of the tubing only
Pump	From pump outlets to pressure transducer inlets <b>(Note:</b> this tubing may already be installed on your workstation)	2 x 100 cm ± 1 cm	PEEK, Orange OD = 1/16" ID = 0.020"	1/4-28 black Delrin fitting on one end of the tubing only
#1	From pressure transducer outlet to dynamic mixer inlet	1 x 28 cm ± 1 cm	PEEK, Orange OD = 1/16" ID = 0.020"	1/4-28 black Delrin fitting on one end of the tubing only
#2	From dynamic mixer outlet to port 5 of AVR7-3 valve	1 x 28 cm ± 1 cm	PEEK, Orange OD = 1/16" ID = 0.020"	1/4-28 black Delrin fitting on one end of the tubing only
#3	From port #4 of AVR7-3 valve to column inlet	1 x 24 cm ± 1 cm	PEEK, Orange OD = 1/16" ID = 0.020"	1/4-28 black Delrin fitting on one end of the tubing only
#4	From column outlet to UV detector inlet	1 x 24 cm ± 1 cm	PEEK, Orange OD = 1/16" ID = 0.020"	1/4-28 black Delrin fitting on one end of the tubing only
#5	From UV detector outlet to conductivity monitor inlet	1 x 15 cm ± 1 cm	PEEK, Orange OD = 1/16" ID = 0.020"	1/4-28 black Delrin fitting on one end of the tubing only
#6	From conductivity monitor to the fraction collector	1 x 52 cm ± 1 cm	Tefzel, clear OD = 1/16" ID = 0.020"	1/4-28 black Delrin fitting on one end of the tubing only

PEEK is a trademark of Victrex, PLC. Tefzel and Delrin are trademarks of E.I. du Pont de Nemours and Company.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [www.bio-rad.com](http://www.bio-rad.com) **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 1 877 89 01 **Belgium** 03 710 53 00 **Brazil** 55 11 3065 7550  
**Canada** 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00  
**France** 01 47 95 69 65 **Germany** 49 89 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300  
**Israel** 03 963 6050 **Italy** 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666  
**New Zealand** 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04  
**Singapore** 65 6415 3188 **South Africa** 27 (0) 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 026 674 55 05  
**Taiwan** 886 2 2578 7189 **Thailand** 1800 88 22 88 **United Kingdom** 020 8328 2000