

BioLogic Maximizer™ Tubing Kit

Catalog #760-2002

The BioLogic Maximizer tubing kit comprises premade inlet tubing to simplify initial setup of your system when a BioLogic Maximizer valve system is used with a BioLogic DuoFlow™ chromatography system workstation containing either F10 or F40 pump heads.

Follow the diagram below to complete the installation.

Note: The PEEK interconnect tubing that connects the Maximizer valves to the workstation pump inlets is supplied separately.

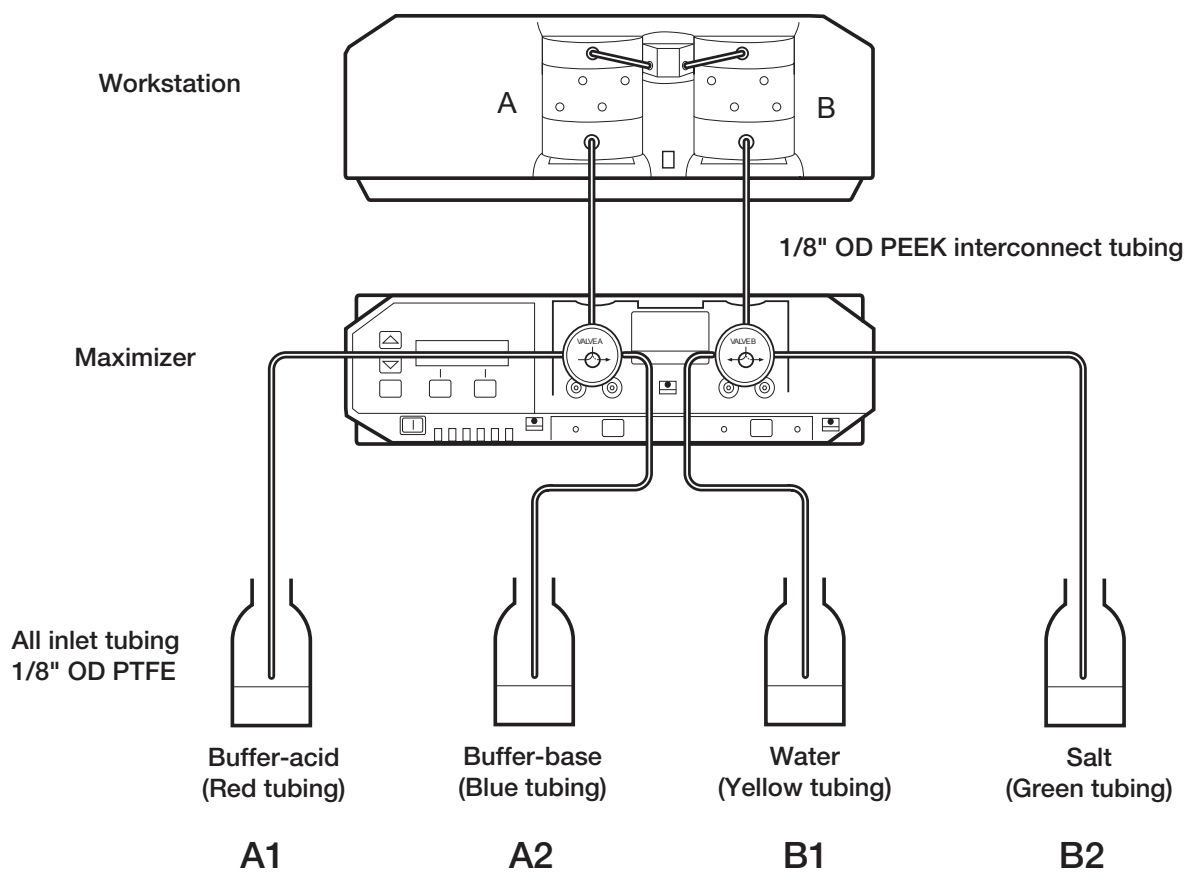


Fig. 1. Plumbing the BioLogic Maximizer valve system with precut tubes from tubing kit.

Table 1. Tubing label and location.

Tubing Label	Tubing Location	Quantity x Tubing Length	Tubing Material, OD/ID	Fittings
Inlet A1	Buffer container to Maximizer valve port A1	1 x 100 cm ± 1 cm	PTFE, red OD = 1/8" ID = 0.062"	1/4-28 black Delrin fitting on one end of the tubing only
Inlet A2	Buffer container to Maximizer valve port A2	1 x 100 cm ± 1 cm	PTFE, blue OD = 1/8" ID = 0.062"	1/4-28 black Delrin fitting on one end of the tubing only
Inlet B1	Water container to Maximizer valve port B1	1 x 100 cm ± 1 cm	PTFE, yellow OD = 1/8" ID = 0.062"	1/4-28 black Delrin fitting on one end of the tubing only
Inlet B2	Salt container to Maximizer valve port B2	1 x 100 cm ± 1 cm	PTFE, green OD = 1/8" ID = 0.062"	1/4-28 black Delrin fitting on one end of the tubing only

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



**Bio-Rad
Laboratories, Inc.**



Life Science
Group

Web site 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