



BioLogic DuoFlow™ pH Monitor

The BioLogic DuoFlow pH Monitor Provides Real-Time
Monitoring of pH Conditions During Sample Runs

BIO-RAD

The BioLogic DuoFlow pH Monitor

The BioLogic DuoFlow inline pH monitor enables real-time monitoring of pH conditions during a sample run, providing you with immediate feedback about the elution profile of your proteins. Key features of the pH monitor include:

- A calomel electrode that, unlike Ag/AgCl electrodes, ensures compatibility with buffers containing Tris
- A single body containing the pH and reference electrodes to minimize reference dry-out and eliminate the need to add electrolyte solutions
- Flow rate capacity of 80 ml/min (75 psi backpressure) to accommodate larger columns
- Biocompatible PEEK flow cell with a swept volume of 80 μ l for high precision and accuracy
- An easy upgrade for any BioLogic DuoFlow system — The pH monitor is supplied with 1/4–28 fitted PEEK tubing for fast and easy addition to any BioLogic DuoFlow system. The electrode connects directly to the BioLogic Maximizer™ valve system's BNC port or, for systems without a Maximizer valve system, via a signal import module (SIM). In addition to being available separately, the pH monitor is included with all BioLogic Maximizer kits, BioLogic DuoFlow Maximizer™ systems, and BioLogic DuoFlow Pathfinder™ systems.



Ordering Information

Catalog #	Description
760-2040	BioLogic DuoFlow pH Monitor , includes SIM module, pH electrode, flow cell, and tubing

Replacement Parts and Accessories

760-2042	pH Electrode
760-2044	pH Flow Cell
760-2046	pH Tubing Kit , includes orange and green PEEK 1/4–28 prefitted tubing lengths for connecting the pH flow cell to the chromatography system
750-0502	Signal Import Module (SIM) , includes 4' communication cable (system cable 17)
760-2034	Universal AC/DC Inline Adaptor for USB Bitbus Device , required when connecting pH monitor or other external detector through SIM module on BioLogic DuoFlow systems

PEEK is a trademark of Victrex PLC.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 424 6723 Australia 61 2 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 31 3689 6600
Canada 905 364 3435 China 86 20 8732 2339 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65
Germany 089 31 884 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300 Israel 03 963 6050
Italy 39 02 216091 Japan 03 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666 New Zealand 0508 805 500
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 27 861 246 723
Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189 United Kingdom 020 8328 2000