



BioLogic DuoFlow System Specifications

Solvent/Buffer Handling

- Dual pump, dual piston, positive displacement solvent delivery module
- 1,000 psi from 0.1–10 ml/min
- High pressure mixing, mixer volume 365 μ l
- Built-in pressure transducer, 34 μ l volume
- Flow accuracy: 0.5% at 1.0 ml/min
2% at 5.0 ml/min
- Flow reproducibility: 0.003 rsd
- Gradient increments: 1%
- Composition accuracy: +0.5% of 100% (>1.0 ml/min)

Detection

- Fixed wavelength, 280/254 nm with auto zero; 0.001–2 AUFS detection range
- 5 mm pathlength analytical flow cell (10 μ l total volume, 6 μ l illuminated volume)
- Optional 2 mm pathlength preparative flow cell (30 μ l)
- Optional 214 nm detection kit (includes lamp, housing and 214 nm filter)
- Optional wavelength filters include: 214, 313, 365, 405, 436, and 546 nm
- Conductivity detection (1 μ S–10 mS; 8 μ l flow cell; accuracy \pm 2% full scale)

Sample Loading

- Fixed sample loops, 25 μ l–5 ml, manual or automated injection
- Large volume loading through the gradient pump. Optional sample inlet valve (SV5-4) available for multiple sample use.
- Software control of peristaltic pump for multiple sample loop fills and large volume load loops

Column Handling

- Single or multiple columns; column switching with optional stream select valves (V9-8 and AV9-8 valves)

Controller

- Internal 486/66MHz computer, 540 Mb, 8Mb RAM hard drive, 3.5" floppy disk
- 14" Super VGA color monitor
- Laboratory-friendly "qwerty" keyboard and mouse

General

- System fittings: 1/4–28 fingertight nuts and ferrules
- System tubing: 1/16" Tefzel (0.062" OD x 0.020" ID)
- Gradient pump inlet tubing: 1/8" OD PTFE
- Dimensions (without collector and recorder): ~68 x 41 x 81 cm (W x D x H), depending upon configuration
- Weight (controller, workstation, monitor, rack, injection valve and detection modules): 34.5 kg/ 76 lbs.
- Wetted materials: 100% biocompatible (PEEK, Tefzel, ceramic)
- Solvent compatibilities: 1 M NaOH, HCl and organic acids, 0.1 M sulfuric acid, 7M urea and GuHCl, 100% MeOH, IPA and acetonitrile, 0.1% TFA, 1% detergents (including SDS, Triton X-100)

Tefzel is a trademark of E.I. du Pont de Nemours and Co.