

## CHROMATOGRAPHY

# C-96 Autosampler

- Easy to connect to the BioLogic DuoFlow chromatography system
- Simple programming via front panel user interface
- Automated, highly reproducible injection of sample volumes from 5  $\mu$ l–5 ml
- Versatile sample handling with three injection modes plus programmable sample and reagent mixing
- Optional sample cooling

## Increase the Efficiency of Your Workflow with the C-96 Autosampler for the BioLogic DuoFlow™ Chromatography System

The C-96 Autosampler enhances the BioLogic DuoFlow chromatography system, providing automated, accurate, and reproducible sample injections. Easy-to-install accessories allow for a wide range of injection volumes from 5  $\mu$ l to 5 ml.

The C-96 Autosampler is available with or without Peltier cooling. The cooling option will maintain heat-sensitive samples at low temperatures during sample handling.

Three injection modes in addition to programmable sample and reagent mixing make the C-96 Autosampler a versatile option for automated sample handling.

### Automated Sample Handling

The C-96 Autosampler handles up to 96 samples. Rapid setup and easy programming add convenient automation to the BioLogic DuoFlow system.

### Biocompatibility

The PEEK fluid path of the C-96 Autosampler ensures biocompatibility for the most sensitive biomolecules.

### Easy Programming

The C-96 Autosampler is fully programmable from a front panel user interface, eliminating the need to learn new software to manage instrument settings. You can program up to nine methods and perform routine maintenance with the touch of a key.



### Easy-to-Install Accessories Expand the Versatility of the C-96 Autosampler

Three dispenser syringes, three sample trays, and a variety of sample loops allow the C-96 Autosampler to accurately and reproducibly perform injections over a wide range of volumes. The optional preparative kit enables injection of up to 5 ml of sample.

### Sample Handling with Minimal Contamination

To prevent contamination between sample injections, the C-96 Autosampler has been designed with buffer tubing between the syringe and injection valves. Washing removes any residual sample from the buffer tubing and sample needles between injections.

**BIO-RAD**

## Specifications

### Instrument

Sample capacity	84 vials of 1.5 ml + 3 reagent vials, 10 ml, for mixing 96 vials of 1.5 ml 24 vials of 10 ml
Vial dimensions (cap included)	Maximum vial height: 48 mm Minimum vial height: 32 mm
Loop volume	50 µl–10 ml
Syringe volume	250 µl, 1,000 µl, 2,500 µl
Optional built-in cooling	4–15°C

### Sampling

Injection modes	Partial loop
	Full loop
	µl pick up
Injection volumes	5 µl–5 ml
Injections per vial	1–9
Programmable methods	9

### Control

BioLogic DuoFlow system	Programmed methods on the C-96 Autosampler are initiated via the BioLogic DuoFlow software
Other systems	Basic control is available via TTL connections

## Ordering Information

Catalog #	Description
760-5010	<b>C-96 Autosampler with Cooling</b> , 110–240 V, includes standard 84+3 vial tray (1.5 and 10 ml), control cable set to connect with BioLogic DuoFlow system, 1 ml syringe, 2 ml sample loop; also includes catalog #760-5014, 760-5026, and 760-0604
760-5011	<b>C-96 Autosampler</b> , 110–240 V, includes standard 84+3 vial tray (1.5 and 10 ml), control cable set to connect with BioLogic DuoFlow system, 1 ml syringe, 2 ml sample loop; also includes catalog #760-5014, 760-5026, and 760-0604
760-5012	<b>Prep Bio Kit</b> , contains 24-position tray for 10 ml vials (22 mm OD), 2.5 ml syringe, 10 ml PEEK loop, 0.75 mm ID PEEK injection valve, prep needle, 6 mm fitting wrench, for use with C-96 autosamplers
760-5013	<b>Syringe</b> , pkg of 1, 1 ml syringe, for use with C-96 autosamplers
760-5014	<b>Connector Kit</b> , contains nuts and ferrules to plumb the syringe valve, for use with C-96 autosamplers
760-5024	<b>Sample Tray</b> , pkg of 1, large-capacity 96-position tray for 1.5 ml vials (12 mm OD), for use with C-96 autosamplers
760-0604	<b>PEEK Tubing</b> , pkg of 1, 1/16" OD x 0.020" ID x 30' high-pressure tubing, rated to 5,000 psi, orange
760-5026	<b>Fittings Kit</b> , contains 10-32 short nuts and ferrules to plumb the injection valve, for use with C-96 autosamplers
760-5027	<b>Needle</b> , pkg of 1, 45 µl needle, for use with C-96 autosamplers
760-5028	<b>Prep Kit Needle</b> , pkg of 1, needle for large sample volume, for use with C-96 autosamplers
760-5017	<b>Analytical Bio Kit</b> , includes 250 µl syringe, 500 µl buffer tubing, 100 and 200 µl PEEK sample loops, standard sample needle, 6 mm fitting wrench, for use with C-96 autosamplers

PEEK is a trademark of Victrex PLC.



**BIO-RAD**

**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 01 877 89 01 Belgium 09 385 55 11 Brazil 55 31 3689 6600  
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 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 Malaysia 60 3 2117 5260 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 3170 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 Thailand 66 2 6518311 United Kingdom 020 8328 2000