



BioFrac™ Fraction Collector

From Discovery to Scale-Up, the BioFrac Fraction Collector
Delivers Flexibility for All Your Application Demands



Reliable and Flexible

Greater Efficiency through a Variety of Collection Racks, Hardware, and Integration



Seamless integration with the NGC™ chromatography system and full compatibility with any chromatography system make the BioFrac fraction collector the full-featured fraction collector for all your separation needs. The BioFrac collector includes two indexed racks, each capable of holding up to 90 tubes (13 mm ID), a diverter valve, and preassembled tubing and extra fittings for both high- and low-pressure systems — everything you need to start collecting immediately!



- 1 Drop Arm** — from deep-well microplates to test tubes, the adjustable drop arm accommodates collection vessel heights up to 150 mm
- 2 Collection Path** — collection flow rates as high as 100 ml/min for preparative applications are time-efficient and practical for processing large-volume samples
- 3 Large Panel Display** — the user-friendly large panel display allows quick programming when building methods; collection options include Collect All, Peak Detection, Time Windows or Volume Windows (up to 20), or a combination of Peak Detection and Time Windows
- 4 Diverter Valve** — included with the BioFrac fraction collector, this valve allows flow diversion and minimizes spillage during tube changes
- 5 Off-the-Shelf Racks** — packaged flat and easily assembled, they can be stored in a drawer to free up precious laboratory space
- 6 Function Keys** — on-the-fly editing allows immediate changes during a method run

A Variety of Collection Racks

The BioFrac collector has the capability to deposit fractions to any of several new custom collection racks, as well as off-the-shelf Scienceware collection racks. Off-the-shelf racks are economical and space-saving and improve workflow by allowing immediate transfer of fractions to cold storage.

Flexible Collection Hardware

BioFrac collector features include reliable serpentine pattern collection and a vertically adjustable drop arm to accommodate various collection vessel heights.

Complete System Integration

The BioFrac collector is easily programmed for simple or complex collection schemes and can be interfaced with a variety of chromatography systems. A 15-pin to bare wire cable is available for non-Bio-Rad pumps, UV detectors, and chart recorders, giving researchers even more flexibility when it comes to application setup.

Versatile Rack Options

The Convenience of Off-the-Shelf Racks

Bio-Rad offers a wide variety of off-the-shelf racking options for collection into test tubes, microtubes, and scintillation vials. Additional Scienceware racks for tubes of various sizes can easily be obtained from providers such as Fisher Scientific and VWR. Scienceware racks are easily assembled and flattened for convenient storage.



The Flexibility of Custom Racks

The optional ice bath/microplate rack offers dual functionality. Designed to protect sensitive samples from degradation, the rack base holds ice while the lid securely holds 12–13 mm tubes. Snap-on insert rings hold tubes firmly in place to allow one-step decanting of noncritical fractions.

Alternatively, the rack lid is designed to securely hold four microplates or racked microtiter tubes when the collection of microliter volumes is required. The accommodation of a variety of microplate well formats offers the flexibility needed for today's researchers.

With the addition of the optional Prep-20 adaptor, up to 20 fractions can be collected to any size bottle, with flow rates up to 100 ml/min for rapid processing.



BioFrac fraction collector with ice bath rack



BioFrac fraction collector with Prep-20 adaptor



The ice bath rack holds up to 4 microtiter trays or microplates

Specifications

Fractionation	Time: 0.01–9,999 min Drop: 1–99,999 drops, flow rate ≤5.0 ml/min Volume: 0.02–99,999 ml
Collection mode	Time, drop, time windows (up to 20 windows), peak detection (threshold), time windows plus peak detection, volume (flow rate ≤100 ml/min)
Peak detection	Threshold, including detection of maxima/minima for collection of double peaks above a set threshold
Detector input	±0–100 mV, ±0–1 V
Wetted parts	All nonmetal, including Tefzel, PEEK
Weight	6.9 kg (15.3 lb)
Dimensions	44.5 x 35.6 x 38.7 cm (17.5 x 14.0 x 15.2")
Test tube change time	~330 ms with 13 mm tubes; ~450 ms with 18 mm tubes
Safety certification	IEC 61010 and CSA 22.2

Ordering Information

Catalog # Description

BioLogic Fraction Collector

741-0002	BioFrac Fraction Collector , 110 V includes power cord, rack set F1, diverter valve, fittings kit, instructions
741-0010	Rack Set F1 (2 x flatpack) , each holds 90 tubes, 12–13 mm diameter, for total of 180 tubes per collection
741-0011	Rack Set F2 (2 x flatpack) , each holds 60 tubes, 15–16 mm diameter, for total of 120 tubes per collection
741-0012	Rack Set F3 (2 x flatpack) , each holds 40 tubes, 18–20 mm, for total of 80 tubes per collection
741-0013	Rack Set H1 (4 x flatpack) , each holds 42 capless 1.5 ml microtubes, for total of 168 microtubes per collection
741-0014	Rack Set H2 (4 x flatpack) , each holds 63 capless 0.5 ml microtubes, for total of 252 microtubes per collection
741-0015	Rack Set H3 (4 x flatpack) , each holds 30 reduced-volume scintillation vials, 16 mm diameter, for total of 120 vials per collection
741-0016	Rack Set H4 (4 x flatpack) , each holds 6 scintillation vials, 30 mm diameter, for total of 24 vials per collection
741-0020	BioFrac H4-High Rack Set (4 x flatpack) , each holds 6 centrifuge tubes, 30 mm diameter, for total of 24 vials per collection
741-0017	BioFrac Ice Bath/Microplate Rack , holds 120 tubes, 12–13 mm diameter; microplate rack feature allows use of up to 4 SBS format microplates in 96-, 48-, 24-, or 12-well configurations; TiterTube® collection can be done using an 8 x 12, 96-tube configuration
741-0018	BioFrac Prep-20 Preparative Rack
741-0007	BioFrac Fraction Collector Fittings Kit , includes replacement fittings and tubing for setup of the fraction collector to the BioLogic™ LP or BioLogic DuoFlow™ chromatography systems
741-0088	BioFrac Microplate Drophead Kit

Tubes

223-9750	Clear Polystyrene Tubes , 13 x 100 mm, 9 ml nominal capacity, 1,000/box
223-9751	Natural Polypropylene Tubes , 13 x 100 mm, 9 ml nominal capacity, 1,000/box
223-9500	Capless Micro Test Tubes , 1.5 ml, polypropylene, natural, for use with micro tube adaptor, 500/box

Cables

731-8263	System Cable 3
731-8286	System Cable 15
731-8287	System Cable 16
731-8290	BioFrac Accessory Cable
731-9009	System Cable 23

PEEK is a trademark of Victrex PLC. Scienceware is a trademark of Bel-Art Products.
Tefzel is a trademark of E.I. du Pont de Nemours and Co.



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 11 5044 5699
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 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** 03 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** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 026674 55 05
Taiwan 886 2 2578 7189 **Thailand** 800 88 22 88 **United Kingdom** 020 8328 2000