

When it comes to
protein purification
hand over
control
to us



The New Econo Gradient Pump Controls
Salt Gradient Formation, Fraction Collection and
Even a Stream-Splitter Valve

BIO-RAD

It's More Than Just a Pump!

The unique Econo Gradient Pump is the one pump you want in your lab to control both your protein purification protocols and new applications based on stream splitting.

For Protein Purification

The Econo Gradient Pump is a 2-channel peristaltic pump perfect for low pressure chromatography applications such as purifying recombinant proteins via affinity resins. For non-affinity tagged proteins, the highly accurate control of isocratic and gradient elution programs means reproducible results, run after run.

FEATURES

- Flow rates of 0.01-20 ml/min per channel, depending on tubing ID
- Control of a Gradient Mixer valve for binary gradient formation
- Simple gradient programming via a 6-step protocol in either Time or Volume units
- One program stored in non-volatile memory
- Purge and Reverse Flow functions
- Control of a diverter valve for one Collection Window
- Stop-start of fraction collectors
- Hold and Pause features available with certain collectors
- Stop-start, pen up-down control of chart recorders
- User-friendly conveniences, including adjustable LCD screen contrast, revolution counter to determine tubing wear
- Uses luer-lock fittings for fast and simple tubing connections
- Compatible with Bio-Rad's BioLogic Duo-Flow chromatography system
- Full range of chromatography media available

To complete your chromatography system, the Econo Gradient Pump may be coupled with Bio-Rad's Model EM-1 UV detector and EG-1 conductivity detector, the Model 1327 chart recorder and either the Model 2110 or Model 2128 fraction collectors (see our Low Pressure Components brochure, bulletin 2437).

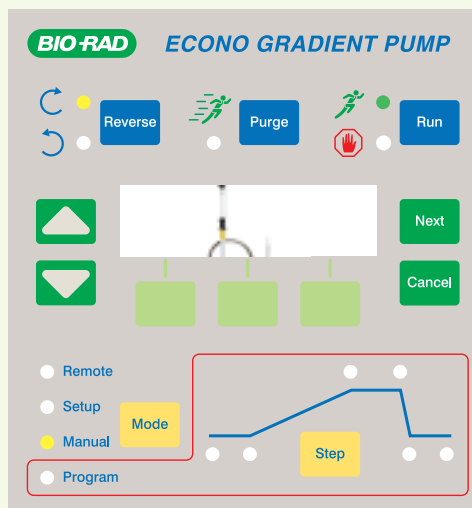
For Stream Splitting Applications

The Econo Gradient Pump controls a new Splitter Valve for stream splitting, on-line bioassays and process sampling applications. It may be used with any chromatography system.

FEATURES

- Splitter Valve with low internal dead-volume of <10 μ l
- Maximum pressure of 30 psi for post-column flow
- With the optional Splitter Valve kit, flow rates as low as 2 μ l/min are achievable
- Variable "% Split" and "Splitter Valve Period" for maximum flexibility and control of your applications

Note that a Diverter Valve and a Splitter Valve cannot be connected to the Econo Gradient Pump at the same time.



Ordering Information

Catalog # Description

ECONO GRADIENT PUMP

- | | |
|----------|--|
| 731-9001 | Econo Gradient Pump, 100/120 V, includes tubing and fittings kit |
| 731-9002 | Econo Gradient Pump, 220 V, includes tubing and fittings kit |

COMBINATIONS

- | | |
|----------|--|
| 731-9030 | Econo Gradient Pump Combo 1, 100/120 V, includes Econo Gradient Pump and Gradient Mixer valve |
| 731-9032 | Econo Gradient Pump Combo 1, 220/240 V, includes Econo Gradient Pump and Gradient Mixer valve |
| 731-9034 | Econo Gradient Pump Combo 2, 100/120 V, includes Econo Gradient Pump, Gradient Mixer valve, MV-6 injection valve and rack with column clamps |
| 731-9036 | Econo Gradient Pump Combo 2, 220/240 V, includes Econo Gradient Pump, Gradient Mixer valve, MV-6 injection valve and rack with column clamps |
| 731-9038 | Econo Gradient Pump Combo 3, 100/120 V, includes Econo Gradient Pump, Gradient Mixer valve and rack with column clamps |
| 731-9040 | Econo Gradient Pump Combo 3, 220/240 V, includes Econo Gradient Pump Gradient Mixer valve and rack with column clamps |

DIVERTER VALVE

- | | |
|----------|----------------------|
| 731-8367 | SVF-3 Diverter Valve |
|----------|----------------------|

SPLITTER VALVE

- | | |
|----------|---|
| 731-9011 | Splitter Valve Kit, includes valve, tubing, adapters, "T" union, back-pressure device |
|----------|---|

CABLES

- | | |
|----------|--|
| 731-9009 | Cable #23, Y-cable connecting Econo Gradient Pump to Model 2128 Fraction Collector and Model 1327 Chart Recorder |
| 731-9010 | Cable #22, Y-cable connecting Econo Gradient Pump to Model 2110 Fraction Collector and Model 1327 Chart Recorder |
| 731-8267 | System Cable 7, 8-pin mini DIN to bare wires for connecting Econo Gradient Pump to generic fraction collectors and chart recorders |

BIO-RAD

**Bio-Rad
Laboratories**

Life Science
Group

Web Site www.bio-rad.com **U.S.** (800) 4BIORAD **Australia** 02 9914 2800 **Austria** (01)-877 89 01 **Belgium** 09-385 55 11 **Canada** (905) 712-2771
China 86-10-62051850/51 **Denmark** 45 39 17 99 47 **Finland** 358 (0)9 804 2200 **France** 01 43 90 46 90 **Germany** 089 318 84-0
Hong Kong 852-2789-3300 **India** (91-11) 461-0103 **Israel** 03 951 4127 **Italy** 39-02-216091 **Japan** 03-5811-6270 **Korea** 82-2-3473-4460
Latin America 305-894-5950 **Mexico** 514-2210 **The Netherlands** 0318-540666 **New Zealand** 64-9-4152280 **Norway** 22-74-18-70 **Russia** 7 095 979 98 00
Singapore 65-2729877 **Spain** 34-91-661-7085 **Sweden** 46 (0)8-55 51 27 00 **Switzerland** 01-809 55 55 **United Kingdom** 0800-181134