



PROCESS-SCALE CHROMATOGRAPHY

Liquid Pump for Inflatable Seals

Liquid Pump for Bio-Rad Process-Scale Chromatography Column Seals

The liquid pump inflates the seals on Bio-Rad process-scale chromatography columns without the need for compressed air. The liquid pump is mounted on top of the column and allows the seal to be manually inflated.

Advantages of the liquid pump:

- Self-contained system eliminates the need for compressed air
- Column stays watertight without permanent air connection
- Manometer is standard equipment

Glycerol is the standard filling for the liquid pump and inflatable seal, but another liquid can be used if specified by the customer.

Specifications

Body and piston	Stainless steel 316L
Seals	EPDM

Dimensions

Liquid pump volume (L)	0.5	0.5	0.5	0.5	0.5	0.5	1.0	1.0	1.0
Column diameter (mm)	100	200	296	350	400	446	596	800	1,000



Ordering Information

Catalog # Description*

Liquid Pump Only

736-0419 **Liquid Pump**, 0.5 L volume
736-0420 **Liquid Pump**, 1 L volume

Column Upgrade with Liquid Pump

736-0981 **Upgrade with Liquid Pump**, 0.5 L volume
736-0982 **Upgrade with Liquid Pump**, 1.0 L volume
736-0983 **Upgrade with Liquid Pump**, 2.0 L volume

* Each catalog number includes the parts necessary for removing the standard air regulator and for mounting and connecting the liquid pump.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 4BIORAD **Australia** 61 02 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 21 3237 9400
Canada 905 364 3435 **China** 86 21 6426 0808 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65
Germany 089 318 84 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 455 8800 **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