



PROCESS-SCALE CHROMATOGRAPHY

Pressure Relief Valve

Protect Columns With Pressure Relief Valve Accessory

The pressure relief valve protects Bio-Rad process-scale chromatography columns and the Bio-Rad® bubble trap from damage during loading of media and column operation. An option for any Bio-Rad process chromatography column, the valve can be attached easily with clamp connections to the top of the column (T-connection) or to the cover of the bubble trap.

Specifications

Valves	Stainless steel 316L, $R_a < 0.4 \mu\text{m}$									
Seals	EPDM									
Clamp connections	Tri-Clamp 25 or 50 (specified by user)									
For column diameter (mm)	100	200	296	350	400	446	596	800	1,000	
Process tube diameter (in)	1/4	3/8	1/2	1/2	3/4	3/4	3/4	1	1	
Clamp connections (mm)	25	25	25	25	25	25	25	50.5	50.5	
For bubble trap	See ordering information (on reverse) to select correct size relief valve based on bubble trap volume.									



Ordering Information

Catalog #	Description*
739-2590	Pressure Relief Valve , for column diameter 100 mm
739-2591	Pressure Relief Valve , for column diameter 200 mm
739-2592	Pressure Relief Valve , for column diameters 296 mm, 350 mm
739-2593	Pressure Relief Valve , for column diameters 400 mm, 446 mm, 596 mm
739-2594	Pressure Relief Valve , for column diameters 800 mm, 1,000 mm
739-2595	Pressure Relief Valve , for bubble trap sizes 0.8 L, 3.5 L, and 5.4 L
739-2596	Pressure Relief Valve , for bubble trap size 9.1 L

* Each catalog number includes one pressure relief valve, one T-connection (for column only), three clamp collars, and three EPDM clamp gaskets.

Tri-Clamp is a trademark of The Ladish Company, Tri-Clover Division.



**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