

PROCESS-SCALE CHROMATOGRAPHY

Pressure Gauge

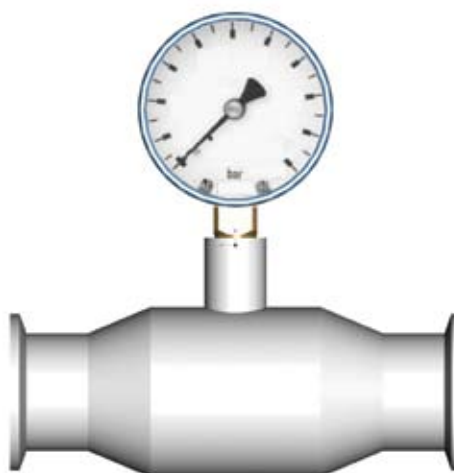
Monitor Column Pressure Easily With Column Accessory

The pressure gauge is used to monitor column pressure on Bio-Rad process-scale chromatography columns. It is equipped with an in-line hygienic separator that does not create a dead leg in the process piping. The pressure gauge is easily attached to the top of the column with clamp connections. Both bar and psi can be read on the gauge indicator.

Specifications

Material									
Pressure gauge									
Wetted parts (tube and diaphragm)									
Connections									
Certificates									
Fill fluid									

For Process-Scale Column Diameter (mm)	100	200	296 350	400 446 596	800 1,000
Pressure gauge tube diameter (in)	1/2	1/2	1/2	3/4	1
Tube length (mm)	114	114	114	114	114
Gauge diameter (mm)	63	63	63	63	63
Clamp connection (mm)	25	25	25	25	50.5
Range (bar)	0-6	0-6	0-6	0-6	0-6



Ordering Information

Catalog #	Description*
739-2360	Pressure Gauge , for column diameter 100 mm
739-2361	Pressure Gauge , for column diameter 200 mm
739-2362	Pressure Gauge , for column diameters 296 mm, 350 mm
739-2363	Pressure Gauge , for column diameters 400 mm, 446 mm, 596 mm
739-2364	Pressure Gauge , for column diameters 800 mm, 1,000 mm

* Each catalog number includes one pressure gauge, two adaptors, two clamp collars, and four clamp gaskets for columns 100 and 200 mm in diameter, and one pressure gauge, two clamp collars, and two clamp gaskets for columns 300 to 1,000 mm in diameter.

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