

PROCESS-SCALE CHROMATOGRAPHY

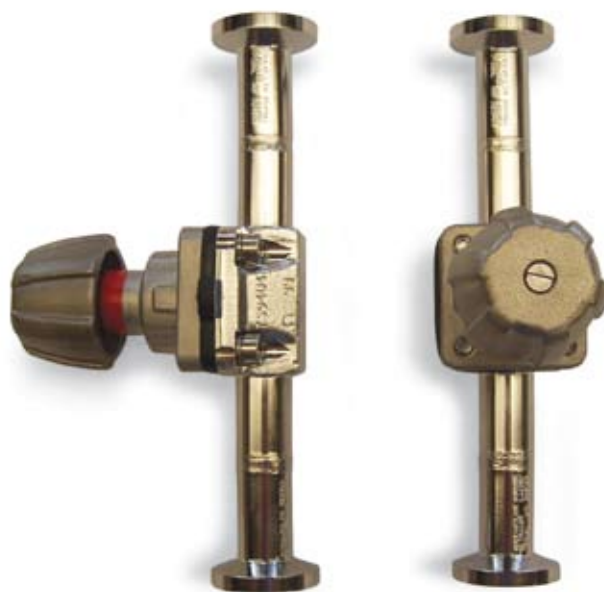
Isolation Two-Way Valves

Take Columns Off-Line With Column Valve Accessory

Isolation two-way valves are attached to the process chromatography column inlet and outlet to isolate the column from the environment during downtime or storage. Isolation two-way valves can also be used for priming the central screw. The valves are attached easily with clamp connections at the top and bottom of the column.

Specifications

Valves	Stainless steel 316L, $R_a < 0.4 \mu\text{m}$								
Valve diaphragms	EPDM								
Clamp connections	Tri-Clamp 25 or 50								
Column diameter (mm)	100	200	296	350	400	446	596	800	1,000
Tube diameter (in)	¼	¾	½	½	¾	¾	¾	1	1
Clamp connections (mm)	25	25	25	25	25	25	25	50.5	50.5
Total length (mm)	129	129	129	129	165	165	165	177	177



Ordering Information

Catalog #	Description*
739-2370	Isolation Two-Way Valve , ¼ in tube diameter
739-2371	Isolation Two-Way Valve , ⅜ in tube diameter
739-2372	Isolation Two-Way Valve , ½ in tube diameter
739-2373	Isolation Two-Way Valve , ¾ in tube diameter
739-2374	Isolation Two-Way Valve , 1 in tube diameter

* Each catalog number includes one valve, two clamp collars, and two 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