



## PROCESS-SCALE CHROMATOGRAPHY

# Bio-Rad® Bubble Trap

### Column Accessory Protects Packed Columns From Air

The Bio-Rad bubble trap prevents air bubbles from entering process chromatography columns. The Bio-Rad bubble trap, which is available in 0.4, 1.7, 2.7, and 4.6 liter sizes, can be used with any Bio-Rad process chromatography column. It is easily mounted on a column and is connected to the column with a flexible hose. An additional benefit of the Bio-Rad bubble trap is that it reduces residual pulsation.

The Bio-Rad bubble trap is delivered with the inlet and outlet connections capped.

#### Specifications

|                                      |   |            |            |            |
|--------------------------------------|---|------------|------------|------------|
| Flanges                              | Stainless steel                                 |            |            |            |
| Tube                                 | Acrylic, glass, or stainless steel              |            |            |            |
| Wetted parts                         | Stainless steel 1.4404, $R_a < 0.4 \mu\text{m}$ |            |            |            |
| Clamp connections                    | Tri-Clamp 25 or 50                              |            |            |            |
| Flexible hose length (mm)            | 1,500   |            |            |            |
| Operating temperature                | 3–60°C  |            |            |            |
| Regulatory certificates              | FDA, 3.1b, and CE                               |            |            |            |
| <b>Bubble trap volume (L)</b>        | <b>0.4</b>                                      | <b>1.7</b> | <b>2.7</b> | <b>4.6</b> |
| Recommended maximum flow rate (L/hr) | 220   | 600        | 1,000      | 3,200      |
| Connection diameter                  |   |            |            |            |
| Upper 1 (in.)                        | 3/4   | 3/4        | 3/4        | 1 1/2      |
| Upper 2 (in.)                        | 3/4   | 3/4        | 3/4        | 1 1/2      |
| Bottom inlet (in.)                   | 3/4   | 3/4        | 1          | 1 1/2      |
| Bottom outlet (in.)                  | 1/2   | 3/8        | 1          | 1          |
| Maximum pressure at 30°C*            |   |            |            |            |
| Glass tube                           | 6 bar   | 6 bar      | 6 bar      | 5 bar      |
| Acrylic tube                         | 6 bar   | 6 bar      | 6 bar      | 5 bar      |
|                                      | 87 psi  | 87 psi     | 87 psi     | 72 psi     |
|                                      | 87 psi  | 87 psi     | 87 psi     | 72 psi     |

\* The bubble trap should be protected from overpressure. A pressure relief valve mounted on the upper clamp is recommended. The Bio-Rad 3/4" pressure relief valve (catalog #739-2595) can be used with 0.4, 1.7, 2.7, and 4.6 liter sizes bubble traps. The Bio-Rad 1" pressure relief valve (catalog #739-2596) can be used with the 4.6 liter bubble traps.





### Ordering Information

| Catalog # | Description*   |
|-----------|--|
| 739-2410  | <b>Bio-Rad Bubble Trap</b> , glass tube, 0.4 L, 220 L/hr flow rate     |
| 739-2412  | <b>Bio-Rad Bubble Trap</b> , glass tube, 1.7 L, 600 L/hr flow rate     |
| 739-2414  | <b>Bio-Rad Bubble Trap</b> , glass tube, 2.7 L, 1,000 L/hr flow rate   |
| 739-2416  | <b>Bio-Rad Bubble Trap</b> , glass tube, 4.6 L, 3,200 L/hr flow rate   |
| 739-2420  | <b>Bio-Rad Bubble Trap</b> , acrylic tube, 0.4 L, 220 L/hr flow rate   |
| 739-2422  | <b>Bio-Rad Bubble Trap</b> , acrylic tube, 1.7 L, 600 L/hr flow rate   |
| 739-2424  | <b>Bio-Rad Bubble Trap</b> , acrylic tube, 2.7 L, 1,000 L/hr flow rate |
| 739-2426  | <b>Bio-Rad Bubble Trap</b> , acrylic tube, 4.6 L, 3,200 L/hr flow rate |

\* Recommended maximum flow rate is shown. Each catalog number includes one bubble trap, four clamp collars, four EPDM clamp gaskets, two clamp caps, and two flexible hoses.

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](http://www.bio-rad.com) **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 31 3689 6600  
**Canada** 905 364 3435 **China** 86 20 8732 2339 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65  
**Germany** 089 31 884 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **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