



PROCESS-SCALE CHROMATOGRAPHY

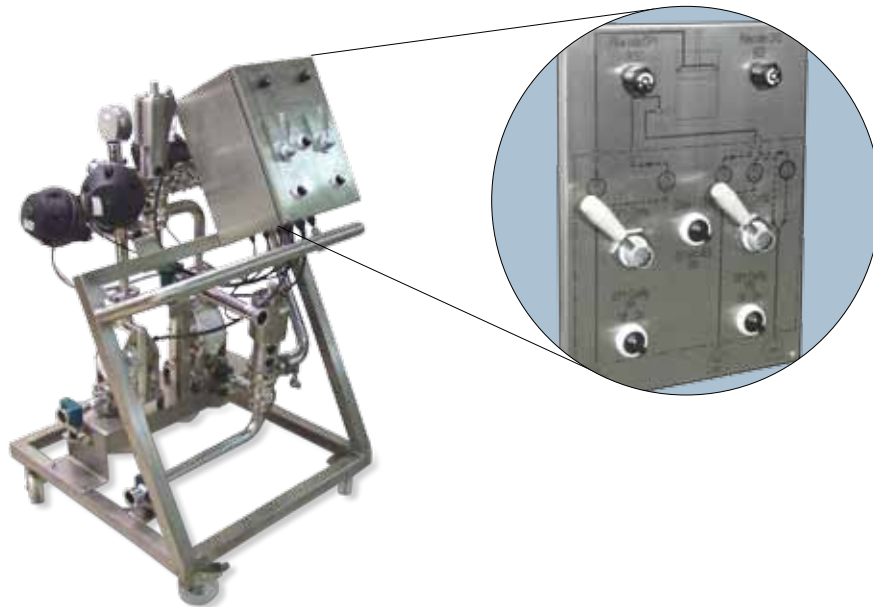
# Bio-Rad® Media Transfer Device

- Low mechanical shear system designed for high value media
- Compact
- Fully pneumatic
- Intuitive user interface
- Complete documentation for regulatory submissions

## The Ideal Counterpart to Bio-Rad® InPlace™ Columns

The Bio-Rad media transfer device is a self-contained movable unit designed to transfer chromatography media between a tank and the column and to deliver cleaning solutions to Bio-Rad InPlace columns for sanitization. The device exploits the full potential of Bio-Rad InPlace column features. Column preparation, unpacking, and cleaning operations are simplified by the device and can be performed in a contained system to preclude contamination.

The Bio-Rad media transfer device is connected to the double manifold of the slurry valves on the Bio-Rad InPlace column and ensures that all wetted surfaces can be sanitized before high-value media are introduced into the column. A dedicated membrane pump gently transfers the media into and out of the column. A second membrane pump is dedicated to transferring the cleaning solutions. The pneumatic control cabinet is semiautomatic. The front panel, comprised of switches and regulators, can be operated intuitively.



An intuitive user interface allows for easy operation of the Bio-Rad media transfer device.



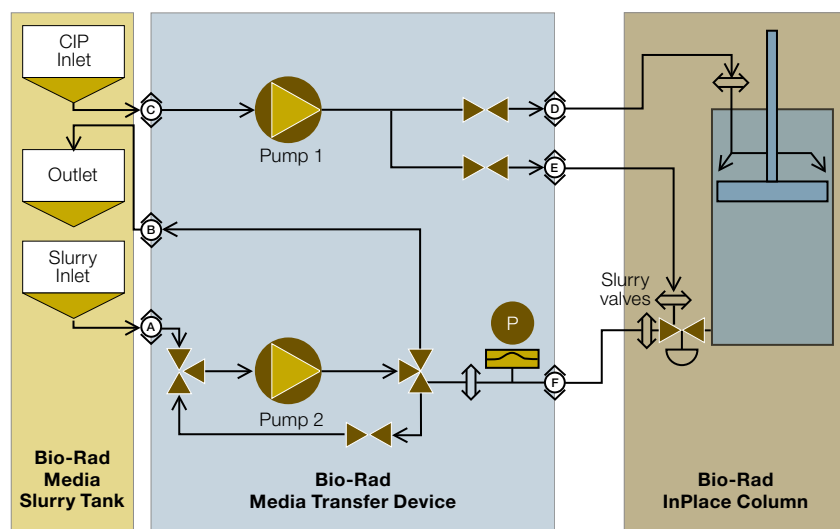
## Manage All Process Operations

With the Bio-Rad media transfer device, you can:

- Flush the slurry valves and manifolds
- Transfer fluid or media from the tank to the column or from the column to the tank
- Clean the column using the integrated cleaning-in-place system of the Bio-Rad InPlace column
- Clean the piping of the device using a clean-in-place protocol

## Regulatory Support and Quality Assurance

The Bio-Rad media transfer device comes with CE certification and is manufactured in our facility, which is certified according to ISO 9001:2008 standards. In Bio-Rad's manufacturing procedures, all steps and components are fully documented. The qualification documentation package assures production in compliance with U.S. and European regulations. The original documentation and one electronic copy are included with each media transfer device. Bio-Rad maintains controlled copies.



Schematic of process setup.

### Specifications

	Flow Rate Range (l/min)		Dimensions (L x W x H), mm	Wetted Surface Finish*
	Min.	Max.		
MTD 50	12	75	740 x 800 x 1,250	$R_a \leq 0.8 \mu\text{m}$
MTD 80	40	150	740 x 800 x 1,250	$R_a \leq 0.8 \mu\text{m}$
MTD 240	118	400	1,000 x 1,100 x 1,700	$R_a \leq 0.8 \mu\text{m}$

\* Higher level of surface finish is available on a special order basis.

### Bio-Rad Media Transfer Device and Column Diameter Compatibility

	Bio-Rad InPlace Column Nominal Internal Diameter, mm						
	400	450	600	800	1,000	1,200	1,400
MTD 50	YES						
MTD 80			YES				
MTD 240						YES	



**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