

PROCESS-SCALE CHROMATOGRAPHY

Bio-Rad® Media Slurry Tank

Optimize Column Operation and Handling of Media

The Bio-Rad media slurry tank is a stand-alone accessory. It can be used with Bio-Rad® EasyPack™ columns combined with a transfer pump, and it can be used with Bio-Rad® InPlace™ columns combined with a Bio-Rad® media transfer device. The tank is available in 100, 300, 500, or 900 L sizes; each has one access lid, one top inlet with clamp connection, and one lower outlet with clamp connection. The tank bottom is sloped for optimized drainability.

The tank uses an agitator with a double propeller to ensure that media are mixed properly and a rotating spray ball for cleaning in place. The tank is designed for use at atmospheric pressure. Material certificates, welding certification, surface finish report, pickling and passivation report, and instruction manual are included.

The Bio-Rad media slurry tank fulfills several functions:

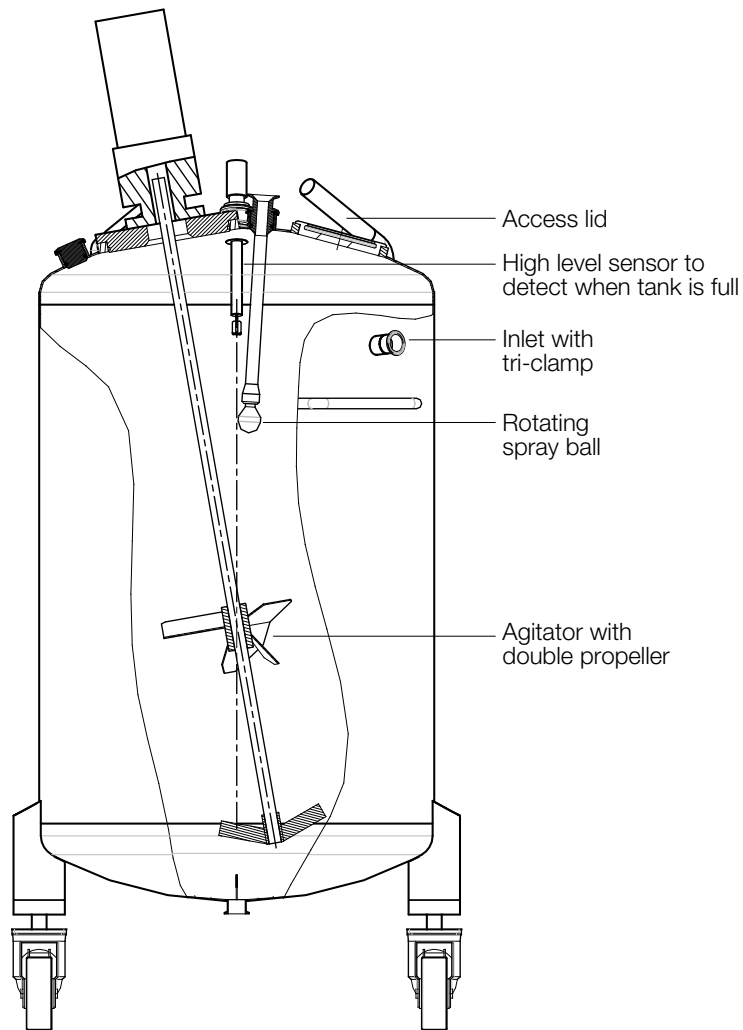
- Resuspends slurry even after complete settlement of the media
- Maintains slurry in suspension during transfer of media to the column
- Rinses interior after media transfer to ensure complete removal of the media
- Cleans and chemically sanitizes tank using spray ball
- Receives resin and rinsing buffer from column during unpacking

The Bio-Rad media slurry tank is also suitable for use with other manufacturers' columns.

Specifications

Tank operating pressure	Atmospheric (<0.5 bar)	
Wetted material		
Parts	Stainless steel 316L	
Seals	EPDM (21 CFR 177.2600)	
Spray ball bushing	PTFE (21 CFR 177.1550)	
Surface finish	R _a ≤ 0.8 μm, mechanically polished; other surface finishes available as options	
Integrated components		
Agitator	Double propeller, 255 r/min maximum	Agitec or equivalent
High level sensor	Vibrating fork	Endress+Hauser or equivalent
Spray ball		Alfa Laval or equivalent
Working pressure	1–3 bar	
Flow rate	80 Lpm maximum	
Variable frequency drive	MICROMASTER 420 or equivalent	Siemens or equivalent





Ordering Information

Catalog #	Description
739-2400	Bio-Rad Media Slurry Tank, 100 L
739-2401	Bio-Rad Media Slurry Tank, 300 L
739-2402	Bio-Rad Media Slurry Tank, 500 L
739-2403	Bio-Rad Media Slurry Tank, 900 L

MICROMASTER is a trademark of Siemens Corporation.



**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