

# Bio-Scale MT Columns



*Adjustable bed support eliminates column voids*

*Easy assembly for packing and cleaning*

*Optimized design of sample applicator for high resolution chromatography*

*Biocompatible materials for preservation of sample biological activity*

*Transparent glass tube allows easy bed inspection*

*Resistance to high pressures allows use of any chromatographic media*

*Fingertight assembly eliminates the need for tools*

The Bio-Scale Columns are available in four sizes (2, 5, 10, and 20 ml) for easy scale up of separation protocols.

## Bio-Scale MT Columns

The Bio-Scale MT columns are suitable for most chromatography applications. The range of columns has been designed specifically to provide high resolution chromatography of biomolecules when packed with any suitable chromatography media including ion-exchange, hydrophobic interaction, hydroxapatite, and affinity supports. The Bio-Scale MT columns provide precise sample application with the low dead volumes required for high resolution separations. The availability of four column sizes (2, 5, 10, and 20 ml) allow easy scale-up of separation and purification protocols.

The Bio-Scale MT column's unique design has been optimized to provide easy packing, bed-height adjustment, sample application and equilibration. The Bio-Scale MT columns offer the chromatographer all the convenience needed when working with the BioLogic System or any medium or high pressure chromatography system.

## Column Specifications

All wetted parts are bio-compatible (Tefzel, polypropylene, polyphenyl sulfide, borosilicate glass and nitrile rubber) and resistant to cleaning and sanitation solvents including 1.0 M NaOH and dilute acetic acid. The column can be used with most common aqueous buffers and certain organics including alcohols and acetonitrile. The Bio-Scale MT columns are not intended for use with halogenated hydrocarbons, aromatic solvents and tetrahydrofuran (THF).

### Ordering Information

Catalog Number	Product Description	Bed Dimension (mm)	Column Volume (ml)	Pressure Limit (psi)
751-0081	Bio-Scale MT2	7 x 52	1.9–2.3	1,000
751-0083	Bio-Scale MT5	10 x 64	4.6–5.7	750
751-0085	Bio-Scale MT10	12 x 88	9.5–11.3	600
751-0087	Bio-Scale MT20	15 x 113	19.4–21.9	500
751-0091	Bio-Scale 2 Replacement Parts Kit, includes 5 frits, 5 distribution screens, 2 o-rings, 1 frit remover			
751-0093	Bio-Scale 5 Replacement Parts Kit, includes 5 frits, 5 distribution screens, 2 o-rings, 1 frit remover			
751-0095	Bio-Scale 10 Replacement Parts Kit, includes 5 frits, 5 distribution screens, 2 o-rings, 1 frit remover			
751-0097	Bio-Scale 20 Replacement Parts Kit, includes 5 frits, 5 distribution screens, 2 o-rings, 1 frit remover			
751-0099	Bio-Scale Fittings Kit, includes 2 Super Flangeless Nuts (1/4 x 28 threads) and 6 ferrules, 2 Flangeless M6 nuts, 4 ferrules and 2 caps, 2 Fingertight II fittings (10–32 threads)			

Note: Each Bio-Scale MT column is supplied with a Bio-Scale Replacement Parts Kit

See the current Bio-Rad catalog for additional information. Specifications are subject to change without notice.



Bio-Rad  
Laboratories

Life Science  
Group

Website [www.bio-rad.com](http://www.bio-rad.com) Bio-Rad Laboratories Main Office 2000 Alfred Nobel Drive, Hercules, CA 94547, Ph. (510) 747-7000, Fx. (510) 747-5800  
Also in: **Australia** Ph. (02-99 74-2000, Fx. (02-99 74-2000) **Austria** Ph. (71-87789 00, Fx. (71) 876 86 29 **Belgium** Ph. (0-385 55 77, Fx. (0-385 65 54  
**Canada** Ph. (905) 772-2777, Fx. (905) 772-2900 **China** Ph. (86-76) 2046522, Fx. (86-76) 2057876 **Denmark** Ph. (39 77 9947, Fx. (39 27 7698  
**Finland** Ph. (00 804 2200, Fx. (00 804 7100) **France** Ph. (7) 43 90 46 90, Fx. (7) 46 77 24 67 **Germany** Ph. (089 37884-0, Fx. (089 37884-700  
**Hong Kong** Ph. 7993300 Fx. 799 1257 **India** Ph. 91-11-467-0103, Fx. 91-11-467-0765 **Israel** Ph. 03 957 4 127, Fx. 03 957 4 129  
**Italy** Ph. 02-27609 7, Fx. 02-27609 399 **Japan** Ph. 03-5877-6270, Fx. 03-5877-6272 **The Netherlands** Ph. 0373 76-540066, Fx. 0373 76-542276  
**New Zealand** Ph. 09-443 3000, Fx. 09-443 3097 **Singapore** Ph. (65) 272-9877, Fx. (65) 273-4838 **Spain** Ph. (91) 667 70 85, Fx. (91) 667 9698  
**Sweden** Ph. 46 (0) 8 627 50 00, Fx. 46 (0) 8 627 54 00 **Switzerland** Ph. 07-809 55 55, Fx. 07-809 55 00  
**United Kingdom** Ph. (0800) 707 734, Fx. 01442 259778