



PROCESS-SCALE CHROMATOGRAPHY Bio-Rad® Packing Motor

Pack Process Chromatography Columns Easily Using Bio-Rad Packing Motors

All Bio-Rad process chromatography columns can benefit from the addition of a Bio-Rad packing motor. A top motor adjusts the height of the adaptor and packs the column via axial compression. Use of the Bio-Rad packing motor eliminates the need to manually activate the central screw. The variable frequency drive of the motor enables highly reproducible axial compression packing.

Motorized axial compression is strongly recommended for columns of diameters between 200 and 1,000 mm; it is compulsory for columns of diameters 1,000 mm and greater.

The motor is controlled by a variable frequency drive installed in an electrical cabinet. The speed of the top piston can be easily adjusted with the potentiometer.

Specifications

Frame and electrical cabinet	Stainless steel 304L
Available power supply	1 AC 220–240 V, 3 AC 220–240 V, 3 AC 380–480 V
Electrical cabinet dimensions	
For column diameters 200–450 mm	400 x 460 x 1,240 mm
For column diameters 600–1,000 mm	600 x 600 x 1,240 mm

Motor and Electrical Cabinet Catalog Number Guide

Column diameter (mm)	200	300	350	400	450	600	800	1,000
Packing motor								
Top	739-2170	739-2171	739-2171	739-2172	739-2172	739-2173	739-2174	739-2175
Electrical cabinet								
Top motor								
1 AC 220–240 V	739-2100	739-2101	739-2101	739-2102	739-2102	739-2103	739-2103	739-2104
3 AC 220–240 V	739-2110	739-2111	739-2111	739-2112	739-2112	739-2113	739-2113	739-2114
3 AC 380–480 V	N/A	739-2121	739-2121	739-2122	739-2122	739-2123	739-2123	739-2124

Note: Please order the appropriate electrical cabinet for the desired motor. For example, if a 3 AC 220–240 V top motor for a 300 mm column is needed, order motor 739-2171 and electrical cabinet 739-2111.



Ordering Information

Catalog # Description

Bio-Rad Packing Motors (Top)

739-2170	Top Motor, for column diameter 200 mm
739-2171	Top Motor, for column diameters 300 mm, 350 mm
739-2172	Top Motor, for column diameters 400 mm, 450 mm
739-2173	Top Motor, for column diameter 600 mm
739-2174	Top Motor, for column diameter 800 mm
739-2175	Top Motor, for column diameter 1,000 mm

Electrical Cabinets

739-2100	Electrical Cabinet for Top Motor, for column diameter 200 mm, 1 AC, 220–240 V
739-2101	Electrical Cabinet for Top Motor, for column diameters 300 mm, 350 mm, 1 AC, 220–240 V
739-2102	Electrical Cabinet for Top Motor, for column diameters 400 mm, 450 mm, 1 AC, 220–240 V
739-2103	Electrical Cabinet for Top Motor, for column diameters 600 mm, 800 mm, 1 AC, 220–240 V
739-2104	Electrical Cabinet for Top Motor, for column diameter 1,000 mm, 1 AC, 220–240 V
739-2110	Electrical Cabinet for Top Motor, for column diameter 200 mm, 3 AC, 220–240 V
739-2111	Electrical Cabinet for Top Motor, for column diameters 300 mm, 350 mm, 3 AC, 220–240 V
739-2112	Electrical Cabinet for Top Motor, for column diameters 400 mm, 450 mm, 3 AC, 220–240 V
739-2113	Electrical Cabinet for Top Motor, for column diameters 600 mm, 800 mm, 3 AC, 220–240 V
739-2114	Electrical Cabinet for Top Motor, for column diameter 1,000 mm, 3 AC, 220–240 V
739-2121	Electrical Cabinet for Top Motor, for column diameters 300 mm, 350 mm, 3 AC, 380–480 V
739-2122	Electrical Cabinet for Top Motor, for column diameters 400 mm, 450 mm, 3 AC, 380–480 V
739-2123	Electrical Cabinet for Top Motor, for column diameters 600 mm, 800 mm, 3 AC, 380–480 V
739-2124	Electrical Cabinet for Top Motor, for column diameter 1,000 mm, 3 AC, 380–480 V



Bio-Rad® InPlace™ column with Bio-Rad packing motor.

BIO-RAD

**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