



## PROCESS-SCALE CHROMATOGRAPHY

# Bio-Rad® Chromatography Station 00

Automated  
preprogrammed  
methods

Minimum cleanroom  
space required

All-in-one  
system design

Controlled by industrial  
wireless tablet PC

U.S. FDA 21 CFR  
Part 11 regulations-  
compliant software

Designed for multiple  
packing methods

Low mechanical shear  
during media transfer

No media loss  
during packing

Adjustable height  
adaptor for axial  
compression

Hoist-free operation

## Automate Process Operations with a Single Unit: Save Cleanroom Space and Time

Bio-Rad's process chromatography station 00 features a high level of automation. The configuration with the motorized Bio-Rad InPlace™ column automatically runs the complete purification process by using preprogrammed methods, starting from column preparation, filling, packing, purification, re-slurrying, and unpacking to cleaning and postprocessing. Operator variances are eliminated and reproducible results are ensured.

The compact design of the Bio-Rad process chromatography station 00 is ideal for small-scale production of pharmaceuticals using column diameters ranging from 100–300 mm, with flow rates from 5–120 L/hr. Its compact size is especially advantageous where space is limited.

### All-in-One Design

The Bio-Rad process chromatography station 00 is composed of the following:

Bio-Rad process skid 00

Bio-Rad process chromatography column

Prep station including components  
for media transfer and column lifting

Wireless tablet PC

Pneumatic or electric components, as well as additional accessories used for media transfer or for lifting column parts, are integrated in the prep station. This design allows the customer to perform all operations for the chromatography process within a single unit, retaining all features and benefits of the separate equipment.

### Controlled by an Industrial Wireless Tablet PC

All actions on the process chromatography station 00 are programmed and controlled with the tablet PC, which is a part of the standard equipment. The wireless connection enables the customer to operate the column and the skid from outside the cleanroom.



**BIO-RAD**

## User-Friendly Software

The software controls and synchronizes the process skid, column motorization, components for media transfer, and accessories for enabling automatic sequences (that is, slurry transfer combined with packing modes or media re-slurrying combined with unloading).

Best practices and methods of Bio-Rad have been created for various chromatography media. These are preprogrammed as recipes in the software and can be used as is or modified to user-defined procedures. This simplifies the column preparation from slurry transfer to unpacking and ensures reproducibility. It also provides required documentation of each step.

The software is designed according to Good Automated Manufacturing Practice (GAMP) 5 guidelines and consists of four access levels (viewer, operator, supervisor, and administrator). It enables the customer to easily program the packing procedure for the column as well as parameters for running the purification on the process skid. Drop-down menus and pop-up screens intuitively guide the user through the programming. Parameter input can be either via keyboard or touch-screen interface of the tablet PC.

Each option includes the required combination of hardware and software modules.

All standard optional components have been validated on the Bio-Rad platform.

### Create Your Process Chromatography Station 00



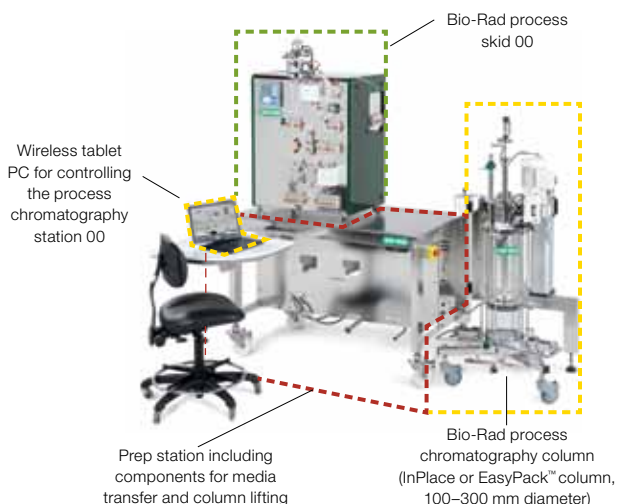
## Ordering Information

Catalog #	Description
<b>Bio-Rad Process Chromatography Skid</b>	
739-3006	Bio-Rad Process Chromatography Skid 00
<b>Select your options for Bio-Rad Process Chromatography Skid 00</b>	
739-3306	3rd pump for sample injection
739-3146	Air sensor after bubble trap
739-3272	Pressure sensor after column
739-3156	Conductivity monitor before column
739-3166	pH monitor before column
739-3126	4 additional outlets

\* Please go to [www.bio-rad.com/process](http://www.bio-rad.com/process) to select the appropriate InPlace or EasyPack column with column IDs between 100 and 300 mm.

For further information, please contact your local Bio-Rad sales representative.

## Components of the Bio-Rad Process Chromatography Station 00



## Modular Concept

The process chromatography station 00 can be adapted to process requirements by integrating options for the Bio-Rad process skid 00 (bulletin 5997) and selecting the appropriate Bio-Rad process column (bulletins 5663 and 5665).

Depending on the column type, the prep station is equipped to drive the following:

Column motorization of the Bio-Rad InPlace or Bio-Rad EasyPack column

Media transfer of the Bio-Rad InPlace column in a contained system

Catalog #	Description
<b>Bio-Rad Process Chromatography Columns</b>	
Varies*	Bio-Rad InPlace Process Column, 180–300 mm
Varies*	Bio-Rad EasyPack Process Column, 100–300 mm
<b>Bio-Rad Prep Station</b>	
739-3810	<b>Configuration 1</b> , InPlace column motorized
739-3820	<b>Configuration 2</b> , EasyPack column motorized
739-3830	<b>Configuration 3</b> , InPlace column nonmotorized
739-3840	<b>Configuration 4</b> , EasyPack column nonmotorized



**BIO-RAD**

**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 21 6169 8500 **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 **Malaysia** 60 3 2117 5260 **Mexico** 52 555 488 7670  
**The Netherlands** 0318 540666 **New Zealand** 64 9 415 2280 **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 3170 **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 **Thailand** 66 2 6518311 **United Kingdom** 020 8328 2000