

CHROMATOGRAPHY

Foresight™ Prepacked Plates and Columns

- Prepacked, ready-to-use formats designed to save process development time
- Different experimental conditions to better define an operational window can be evaluated
- High-throughput experiments with minimal sample requirements can be performed
- Available in a variety of chromatography media modes designed for large-scale bioprocess
- Compatible with robotic liquid handling workstations

Speed and Convenience for Process Developers

Product Information

Foresight plates and columns are prepacked with a range of Bio-Rad's process chromatography media, offering process scientists convenience and reliability for their high-throughput (HT) experimentation needs. Their robust design allows process scientists to use the prepacked formats through the entire purification development cycle, from HT media screening to small-scale methods development and scale-up optimization.

Foresight Plates

Foresight plates are 96-well filter plates prepacked with chromatography media. These plates, with each well prefilled with 20 µl of media, can be used for HT evaluation of various binding, elution, or wash conditions for different chromatography media. Foresight plates can be processed using

manual or automated robotic systems, and sample collection may be achieved through centrifugation or vacuum manifold techniques.

Foresight RoboColumn Units

Foresight RoboColumn units are prepacked microcolumns used for HT process development. Their design is fully compatible with robotic liquid handling systems, allowing process developers to perform automated HT parallel media screening and/or small-scale method development. Foresight RoboColumn units are offered as single rows with each row containing eight columns. They are available in two volumes — 200 µl or 600 µl. Multiple rows up to 12 can be combined to form a 96-column array, allowing users to evaluate various chromatography media and conditions in parallel.

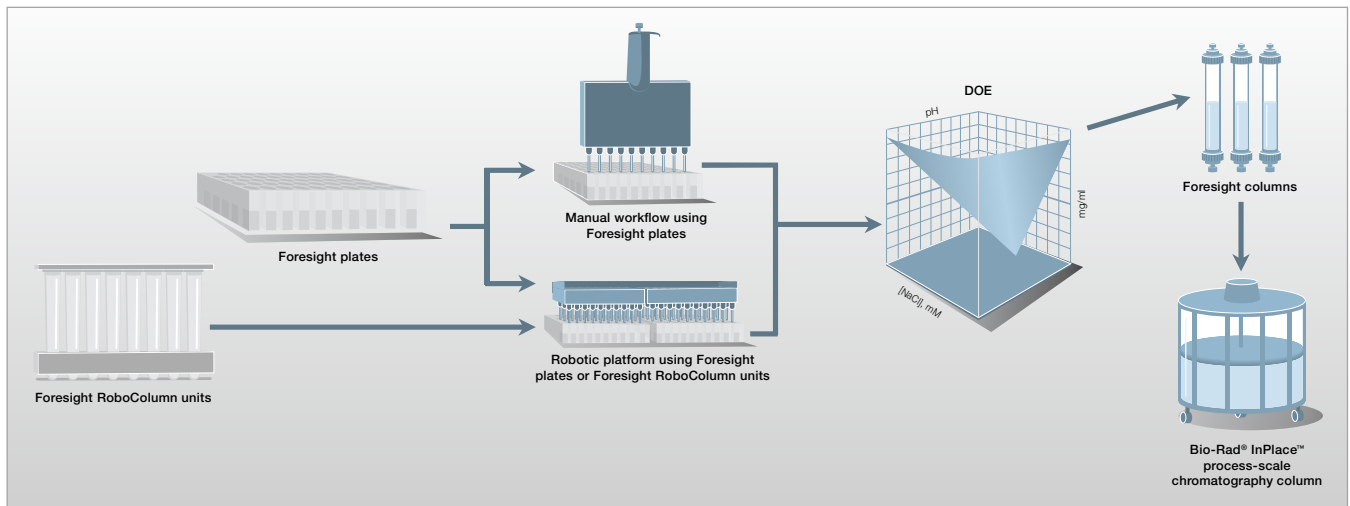


Fig. 1. Purification scale-up workflow shows the use of HT screening and DOE to define an operational window for biomanufacturing.



Foresight Columns

Foresight columns, available in 1-ml or 5-ml formats, may be used for qualitative media screening experiments and/or method development. Foresight columns have standard HPLC connections, which make them compatible with the NGC™ chromatography system and other liquid chromatography systems. These columns are ideal for confirming the operational window obtained through design of experiments (DOE) in trials with Foresight plates and/or Foresight RoboColumn units.

Applications and Regulatory Support

Regulatory support files and application notes are available upon request. Bio-Rad Laboratories is an ISO 9001 registered corporation.

For more information about prepacked columns and plates, please visit www.bio-rad.com/foresight.

Ordering Information

Catalog # Description

Foresight Plates*

732-4714	Foresight™ UNOsphere™ Q Plates, 20 µl
732-4710	Foresight UNOsphere S Plates, 20 µl
732-4712	Foresight UNOsphere Rapid S Plates, 20 µl
732-4703	Foresight™ Nuvia™ Q Plates, 20 µl
732-4701	Foresight Nuvia S Plates, 20 µl
732-4705	Foresight™ Nuvia™ cPrime™ Plates, 20 µl
732-4716	Foresight™ CHT™ Type I Plates, 40 µm, 20 µl
732-4718	Foresight CHT Type II Plates, 40 µm, 20 µl

* Package size: 2 x 96-well plates.

Foresight RoboColumn Units**

732-4819	Foresight UNOsphere Q RoboColumn Units, 200 µl
732-4820	Foresight UNOsphere Q RoboColumn Units, 600 µl
732-4813	Foresight UNOsphere S RoboColumn Units, 200 µl
732-4814	Foresight UNOsphere S RoboColumn Units, 600 µl
732-4816	Foresight UNOsphere Rapid S RoboColumn Units, 200 µl
732-4817	Foresight UNOsphere Rapid S RoboColumn Units, 600 µl
732-4804	Foresight Nuvia Q RoboColumn Units, 200 µl
732-4805	Foresight Nuvia Q RoboColumn Units, 600 µl
732-4801	Foresight Nuvia S RoboColumn Units, 200 µl
732-4802	Foresight Nuvia S RoboColumn Units, 600 µl
732-4807	Foresight Nuvia cPrime RoboColumn Units, 200 µl
732-4808	Foresight Nuvia cPrime RoboColumn Units, 600 µl
732-4822	Foresight CHT Type I RoboColumn Units, 40 µm, 200 µl
732-4823	Foresight CHT Type I RoboColumn Units, 40 µm, 600 µl
732-4825	Foresight CHT Type II RoboColumn Units, 40 µm, 200 µl
732-4826	Foresight CHT Type II RoboColumn Units, 40 µm, 600 µl

** Package size: one row of eight columns.



Foresight Columns

732-4732	Foresight UNOsphere Q Column, 1 ml
732-4752	Foresight UNOsphere Q Column, 5 ml
732-4730	Foresight UNOsphere S Column, 1 ml
732-4750	Foresight UNOsphere S Column, 5 ml
732-4731	Foresight UNOsphere Rapid S Column, 1 ml
732-4751	Foresight UNOsphere Rapid S Column, 5 ml
732-4721	Foresight Nuvia Q Column, 1 ml
732-4741	Foresight Nuvia Q Column, 5 ml
732-4720	Foresight Nuvia S Column, 1 ml
732-4740	Foresight Nuvia S Column, 5 ml
732-4722	Foresight Nuvia cPrime Column, 1 ml
732-4742	Foresight Nuvia cPrime Column, 5 ml
732-4735	Foresight CHT Type I Column, 40 µm, 1 ml
732-4755	Foresight CHT Type I Column, 40 µm, 5 ml
732-4736	Foresight CHT Type II Column, 40 µm, 1 ml
732-4756	Foresight CHT Type II Column, 40 µm, 5 ml

Chromatography Media Available in Foresight Formats

Chromatography Media	Mode
UNOsphere Q	Strong anion
UNOsphere S	Strong cation
UNOsphere Rapid S	Strong cation
Nuvia Q	Strong anion
Nuvia S	Strong cation
Nuvia cPrime	Mixed mode Hydrophobic: cationic
CHT Type I — 40 µm particle size	Mixed mode Metal affinity: cationic
CHT Type II — 40 µm particle size	Mixed mode Metal affinity: cationic

RoboColumn is a trademark of Atoll GmbH.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site 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 11 5044 5699
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 9532 220 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 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 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 Thailand 800 88 22 88 United Kingdom 020 8328 2000