

RoboColumn[®] Formats for HTPD

Prepacked Columns
for Resin Screening and
Method Development



Benefits

- Column format and high quality packing enables mimicking of all steps of a chromatography separation in dynamic mode. Data obtained are comparable to those obtained in larger columns
- Miniaturized format facilitates screening of chromatography resins, reduces sample consumption, and saves time
- Perfectly fitted to Design of Experiment (DoE) requirements, allowing a large amount of experimental work in a minimum of time
- 96-well format allows easy integration with automation systems for High Throughput Process Development (HTPD). Multiple parameters can be tested in parallel, e.g., residence time

Product Information

RoboColumns are miniaturized chromatography columns pre-packed with chromatography resins. They are designed for fully automated and parallel chromatographic separations with robotic liquid handling workstations, such as the Freedom EVO[®] from Tecan.

RoboColumns are packed with ion exchange, mixed-mode, affinity or any other chromatography resins upon request. The column units are packed by Repligen[™] and are identical to OPUS RoboColumns. Two formats are available, 200 μ L and 600 μ L, with a bed height of 10 and 30 mm respectively.

Description

The 200 µL column volume is recommended for resin and process conditions screening, while the 600 µL is used for purification process optimization or when a residence time above 4 minutes is required (e.g., capacity study on mixed-mode resins).

The columns are supplied as rows of eight pre-packed units, with two removable silicon cover seals for proper storage. A 96-well array plate is available to arrange up to 96 RoboColumn units, and the user makes the column selection as needed. The miniturized columns can be reused.



Technical Data

RoboColumns

Column volume	200 µL	600 µL
Bed height	10 mm	30 mm
Column inner diameter	5 mm	
Column material	Polypropylene	
Chemical stability	All commonly used aqueous buffers, pH 1 to 14	
Avoid	Halogenated organic solvents, hexane	
Storage solution	20% ethanol 1 M NaCl	
Recommended storage temperature	2 to 8 °C	
Working temperature	4 to 30 °C	
Maximum working pressure	Up to 8.0 bar g (116 psi g)	
Flow rate	16 to 1,000 cm/hr ⁽¹⁾	
Laser label	Placed on every individual column body, and containing: Product number Resin name Column volume in µL	

96-Well Array Plate

Size	128.3 × 86.0 × 14.0 mm
Material	Polyoxymethylene copolymer
Number of wells	96

(1) The Liquid Transfer menu of the Te-Chrom™ Wizard enables set up of the flow rate for each individual dispensing and chromatography step (5.1.1.6 Liquid Transfer, Te-Chrom Wizard Software Manual, 396076, en, V1.0).

Applications

RoboColumns are suitable for a large variety of applications such as:

- Parallel screening of chromatography resins
- Parallel screening and optimization of chromatographic conditions in downstream process development (conductivity, pH)
- Scale down experimental work
- Determination of dynamic binding capacity
- Determination of the optimal residence time
- Optimization of the regeneration step

Application Example: Development of the Capture Step of a Monoclonal Antibody (MAb) Expressed in a Crude CHO Feedstock

MAb binding conditions were tested on 3 mixed-mode resins and dynamic binding capacity (DBC) was evaluated based on breakthrough (BT) curve (Figure 1).

The RoboColumn was overloaded with MAb CHO feedstock at 0.52 mg/mL and flowthrough (FT) was collected as one-column-volume (CV) fractions (200 μ L each) in 350 μ L 96-well plate. The amount of MAb recovered in the FT fractions was evaluated using Bio-Layer Interferometry (BLI) with protein A sensor (Octet[®] System, ForteBio). It was plotted against the volume of feedstock loaded per RoboColumn unit. 1/10 of the MAb concentration (C/10) was 0.052 mg/mL (52 μ g/mL in Figure 1). The volume equivalent at 10% BT (Veq10%BT) was the volume of MAb loaded when reaching C/10 in FT fractions (8.56 mL in Figure 1).

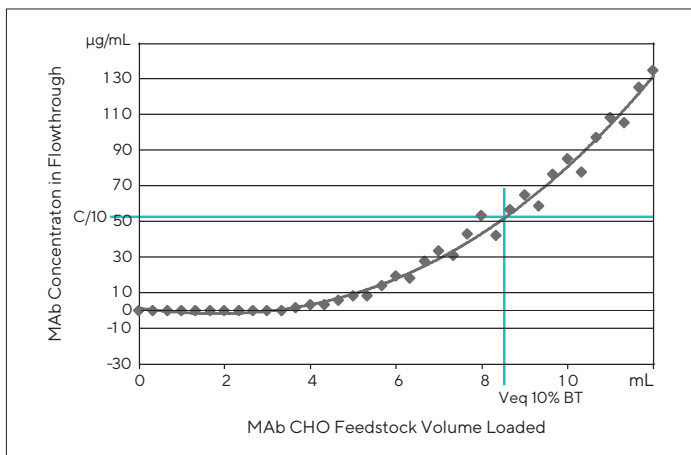


Figure 1: Breakthrough Curve on MEP HyperCel Resin in 200 μ L RoboColumn. Breakthrough during loading of MAb CHO feedstock (0.52 mg/mL) at ~ 4 minute RT.

DBC10% BT (mg/mL resin) =

$$\frac{\text{Veq10\%BT (mL)} \times \text{MAb concentration (mg/mL)}}{\text{Column volume (mL)}}$$

* void volume was taken into account

Using this breakthrough curve method, the DBC for MEP, HEA and PPA HyperCel resins was determined on 200 μ L RoboColumns (Figure 2).

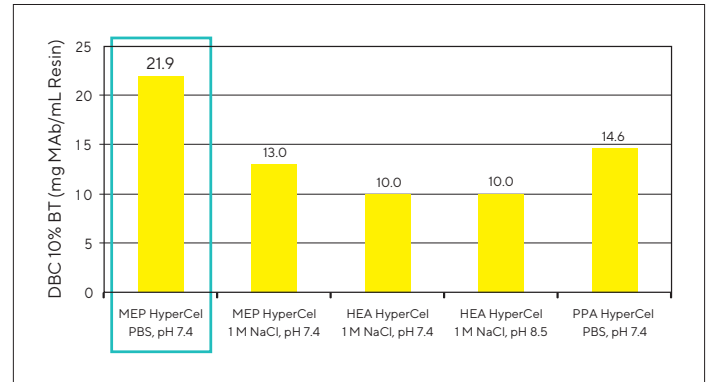


Figure 2: DBC on 3 Mixed-Mode HyperCel Resins in RoboColumns for Different Binding Conditions

This comparison showed that MAb binding was maximized on MEP HyperCel resin in PBS, pH 7.4. The selected condition was then transferred to 1 mL PRC column for confirmation of the DBC, approximately 21 mg/mL (Figure 3).

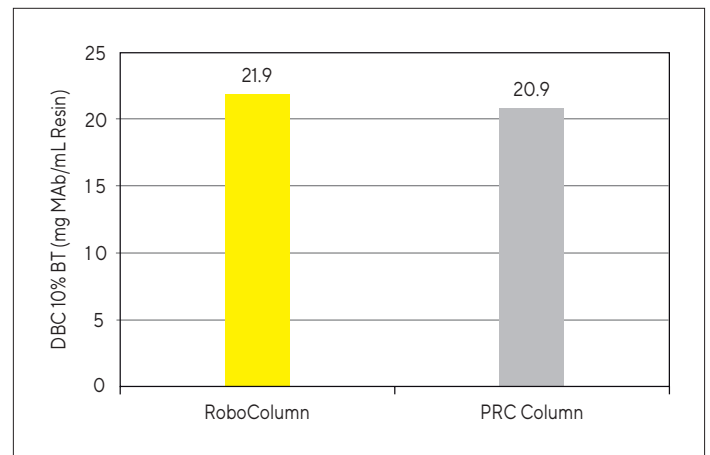


Figure 3: DBC on MEP HyperCel Resin in RoboColumn and in PRC Column

A 335-fold scale-up of the MAb capture step was finally carried out from 200 μ L RoboColumn to a lab-scale column (2.5 cm I.D) packed with 67 mL MEP HyperCel resin (Figure 4).

The analysis of fractions obtained on RoboColumn and a 2.5 cm ID lab-scale column were in the same magnitude. Comparable data in terms of recovery (100%) and purity (90%) were evidenced (Figure 5, Table 2).

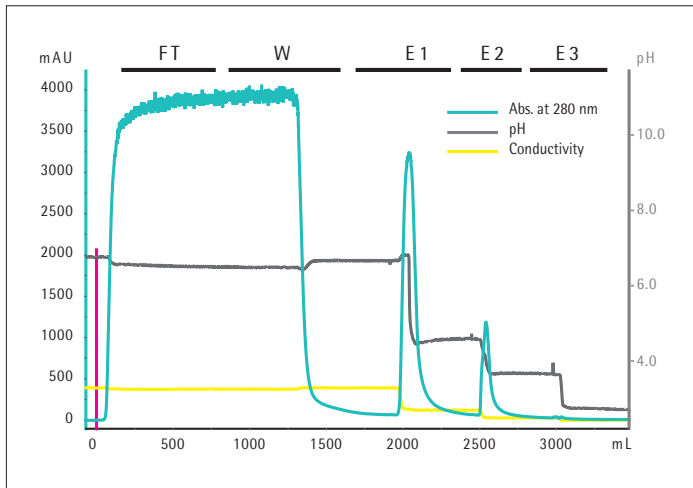


Figure 4: 335-fold Scale-Up of the MAb Capture Step from RoboColumn to a 2.5 cm ID Lab-Scale Column

Chromatogram of MAb Capture Step on a 2.5 cm ID lab-scale column (67 mL). Equilibration in PBS, pH 7.4. Loading of MAb feedstock = 21 mg MAb/mL MEP HyperCel resin. Residence time = 4 minutes; FT: Flowthrough; W: Wash; E1, E2 and E3: Elution at pH 5, 4 and 3 respectively.

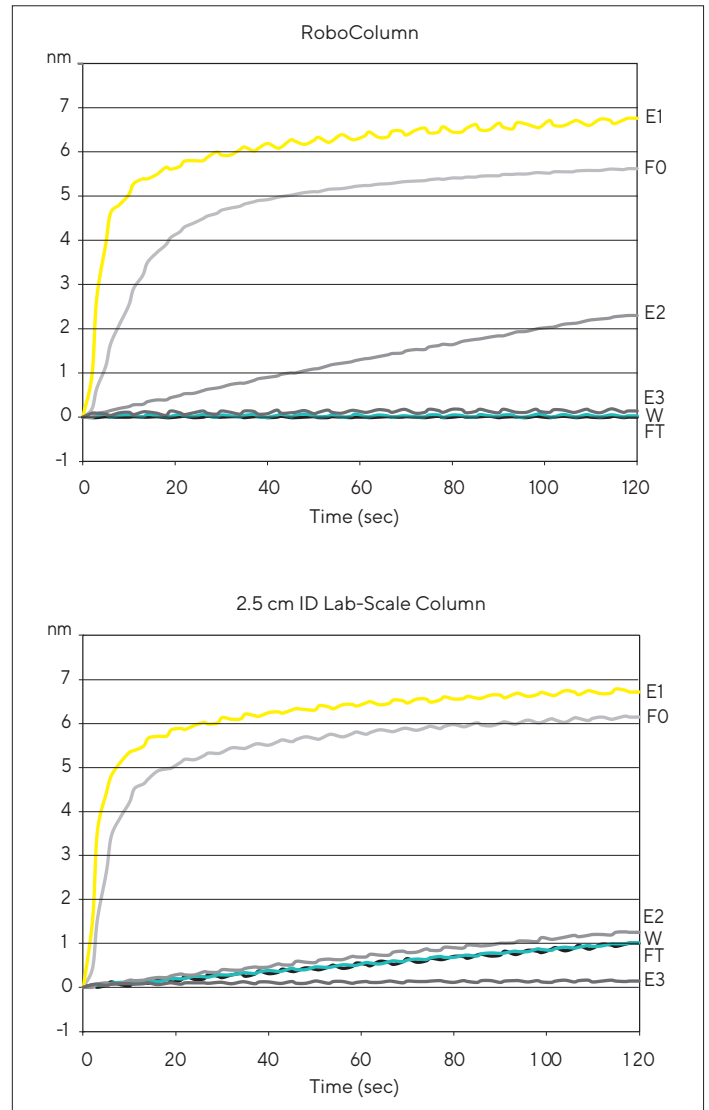


Figure 5: Fraction Analysis From RoboColumn and a 2.5 cm ID Lab-Scale Column Using BLI

Interaction of the MAb with protein A sensor was measured using BLI. Signal is correlated to MAb concentration using a standard curve (data not shown). Residence time = 4 minutes. FO: Sample load; FT: Flowthrough; W: Wash; E1, E2 and E3: Elution at pH 5, 4 and 3 respectively.

Table 2: Summary of the MAb Capture Step on RoboColumn or a 2.5 cm ID Lab-Scale Column

Column Type	Volume (mL)	MAb Loaded (mg)*	Loading (mg MAb/mL resin)	Recovery (%)*	Purity (%)**
RoboColumn	0.2	4	22	102	86
Lab Scale Column 2.5 cm I.D.	66.7	1376	21	99	91

* Evaluated using BLI technology (Octet System, ForteBio)

** Evaluated using LabChip® GXII technology (PerkinElmer)

Handling

RoboColumns are designed for robotic handling. When using the high-throughput process development approach for RoboColumns, it is suggested to use Design of Experiments (DoE) criteria for the experimental set-up. This enables screening of many different chromatographic parameters such as resin type, pH, conductivity| ionic strength, etc., efficiently and simultaneously.

For experimental set-up and the configuration of the 96-well array plate, the Te-Chrom Wizard is recommended. The Te-Chrom Wizard is a dialog-based graphic user interface offered by Tecan and based on the Freedom EVOware software. The Te-Chrom Wizard was developed for the use of RoboColumns and provides the possibility to configure hardware, plate layout, and the chromatographic process as well as to set all process relevant chromatographic parameters, e.g., flow rate, volume without direct script writing.

Ordering Information

Description	Part Number
RoboColumn MEP HyperCel 200 µL, row of 8	SR2MEP
RoboColumn MEP HyperCel 600 µL, row of 8	SR6MEP
RoboColumn HEA HyperCel 200 µL, row of 8	SR2HEA
RoboColumn HEA HyperCel 600 µL, row of 8	SR6HEA
RoboColumn PPA HyperCel 200 µL, row of 8	SR2PPA
RoboColumn PPA HyperCel 600 µL, row of 8	SR6PPA
RoboColumn HyperCel STAR AX 200 µL, row of 8	SR2STARAX
RoboColumn HyperCel STAR AX 600 µL, row of 8	SR6STARAX
RoboColumn CM Ceramic HyperD F 200 µL, row of 8	SR2CMCHDF
RoboColumn CM Ceramic HyperD F 600 µL, row of 8	SR6CMCHDF
RoboColumn CMM HyperCel 200 µL, row of 8	SR2CMM
RoboColumn CMM HyperCel 600 µL, row of 8	SR6CMM
96-well RoboColumn array plate	SR96WAP

¹ Repligen is a trademark of Repligen GmbH


² Te-Chrom™ is a trademark of Tecan

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178

 For further contacts, visit
www.sartorius.com