



Revision history: Pub. No. 100069846 F

Revision	Date	Description
F	14 March 2024	Updated: Resin characteristics Table, Documentation Table, and Ordering Table. Added topic Connecting the column.
E	14 May 2020	Add Oligo dT(25) information. Update manufacturer address.
D	25 February 2019	Table 1: Change chemistry for HS and XS from strong anionexchange to strong cation exchange, change HQ to stronganion exchange, remove HE50. Ordering: Delete HE.
C	6 March 2018	Add affinity chromatography products and 0.8cmD x 10cmL columns.
B	5 March 2018	Change document to small format.
A	31 October 2017	New document. Add HIC products. Replaces 4482551.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2017-2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

For support visit [thermofisher.com/support](https://thermofisher.com/support)  
thermofisher.com  
14 March 2024

POROS™ GoPure™ Pre-Packed Chromatography Columns

Pub. No. 100069846 Rev. F

**WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from [thermofisher.com/support](https://thermofisher.com/support).

Product information

POROS™ GoPure™ pre-packed columns are for screening of POROS™ resins or small scale purification.

Table 1 Resin characteristics

POROS™ Resin	Chemistry	Surface functionality
50 HS	Strong cation exchange	Sulphopropyl
XS		
50 PI	Weak anion exchange	Polyethyleneimine
50 D		Dimethylaminopropyl
50 HQ	Strong anion exchange	Quaternized polyethyleneimine
XQ		Proprietary quaternary amine
Ethyl	Hydrophobic interaction	Novel ethyl
Benzyl		Low-density benzyl/aromatic
Benzyl Ultra		High-density benzyl/aromatic
Caprylate	Mixed-mode cation exchange	Caprylic (octanoic) acid
Oligo dT(25)	Affinity	Poly-deoxythymidine (dT25-mer) ligand
MabCapture™ A		Protein A ligand (outside supplier)
AAVX <sup>[1]</sup>		AAVX ligand
AAV8 <sup>[1]</sup>		AAV8 ligand
AAV9 <sup>[1]</sup>		AAV9 ligand

[1] POROS™ CaptureSelect™

Specifications

Table 2 Column specifications

Condition	Specification
Dimension (nominal volume)	0.5 cmD x 1cmL (0.2 mL), 0.5 cmD x 5 cmL (1 mL), 0.8 cmD x 10 cmL (5 mL)
Maximum pressure	30 bar <ul style="list-style-type: none"><li>Higher pressures will be observed during removal of the ethanol storage solution (1–3 CV)</li><li>POROS™ XQ: Higher column backpressure may be observed with low-conductivity buffers and solutions(&lt;5mS/cm)</li></ul>
Recommended flow rate for process development	≥3 min residence time (0.33 mL/min on a 1-mL column or 1.67 mL/min on a 5-mL column, equates to 400 cm/hr on a 20 cmL column)
Column tube and end cells	High density polyethylene (HDPE)
Frits	Polyethylene/polypropylene 14 µm
Connections	Fingertight for 1/16-inch tubing, female (10-32)
Shipping solution	<ul style="list-style-type: none"><li>AEX and CEX Columns: 20% EtOH, 150 mM NaCl</li><li>HIC and Affinity Columns: 20% EtOH</li></ul>
Asymmetry	0.80 to 1.60
Efficiency	2,500–4,000 plates/m depending on the resin used

For Research Use Only. Not for use in diagnostic procedures.

Guidelines for connecting columns

- Connect the column to the chromatography system using fingertight 1/16", 10-32 fittings.
- Take care not to over-tighten or cross-tread the fittings as this can cause damage to the end fitting.
- When connecting or disconnecting the column, hold the column end-fitting, not the column body to avoid accidentally loosening the fitting.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have any questions, please contact Life Technologies at [www.thermofisher.com/support](http://www.thermofisher.com/support).

Documentation

Publication	Pub. No.
POROS™ Anion Exchange Resins: XQ, 50 HQ, 50 PI and 50 D Product Information Sheet	100025569
POROS™ Strong Cation Exchange Resins: XS and 50 HS Product Information Sheet	100031321
POROS™ 50 HE Product Information Sheet	4327859
POROS™ Ethyl, Benzyl and Benzyl Ultra Hydrophobic Interaction Chromatography Resins Product Information Sheet	100063752
POROS™ Oligo [dT]25 Affinity Resin Product Information Sheet	100092591
POROS™ Protein A Affinity Resins: MabCapture™ A and MabCapture™ A Select User Guide	4383704
POROS™ CaptureSelect™ AAV Resins: AAV8, AAV9, AAVX User Guide	100038399
POROS™ Caprylate Mixed-Mode Chromatography Resin User Guide	MAN0029949

Ordering information

POROS™ GoPure™ Pre-packed column	Cat. No.		
	0.5cmD × 5cmL, 1 mL	0.8cmD × 10cmL, 5 mL	0.5cmD × 1cmL, 0.2 mL
50 D	4481319	A36642	—
50 HS	4481316	A36637	—
50 HQ 50	4481315	A36639	—
MCA	A36644	A36643	—
MCAS	A36646	A36645	—
50 PI	4481318	A36641	—
XQ	A25812	A36640	—
XS	4481317	A36638	—
Benzyl	A34984	A36654	—
Benzyl Ultra	A34985	A36655	—
Ethyl	A34983	A36653	—
AAVX	A36652	A36651	—
AAV8	A36648	A36647	—
AAV9	A36650	A36649	—
Oligo dT[25]	A48352	A48351	A48607
Caprylate	A51054	A51055	—