

Got Purification Troubles?

Improve mAb Purification Cost Efficiency With the Right Resin



The Challenge

Are you trying to improve your process economics? Even after selecting mixed-mode resins with reasonable dynamic binding capacities (DBC), are you unable to lower your process purification costs?



Purification Tip

Only a long-lasting resin with consistent purification performance provides the desired process consistency and economics. Check how many cycles your resin lasts for and if it gives reproducible results at all cycles. The performance of some resins deteriorates over several cycles while others have a short life cycle.



Our Approach

Cycling of a protein mixture (approx. pI 4.5–9.5) and mAb G (approx. pI 9.2–9.5) on two commercially available hydroxyapatite resins showed that Bio-Rad's CHT Ceramic Hydroxyapatite XT Media lasted for over 100 cycles, which was about three times more than the other resin (Table 1). In addition, CHT XT showed the same retention time and peak width even in the 102nd cycle (Figure 1A), whereas a shift in retention time and peak broadening was seen with the other resin in the 27th cycle (Figure 1B). Refer to Bio-Rad bulletin 7080 for technical details on CHT XT.

Table 1. Life of CHT XT versus Ca⁺⁺Pure-HA.

Media	Completed Cycles
CHT XT	102
Ca ⁺⁺ Pure-HA	32

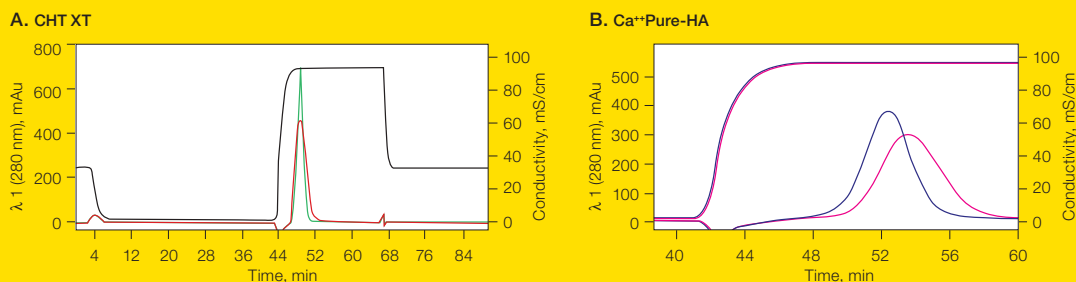


Fig. 1. Purification of mAb G. **A.** on CHT XT there is no change in the retention time or peak width, which leads to reproducible purification. CHT XT cycle 2 (—); CHT XT cycle 102 (—). **B.** on Ca⁺⁺Pure-HA there is a broader peak and delayed elution time at cycle 27. Ca⁺⁺Pure-HA cycle 2 (—); Ca⁺⁺Pure-HA cycle 27 (—).

Go to bio-rad.com/CHTXT and explore the various features of CHT XT.



BIO-RAD

Innovative Products. Customized Solutions.

CHT Ceramic Hydroxyapatite XT Media

CHT XT will put you on your path to:

- Efficient, single-step clearance of impurities
- Great process economics (cost and time efficiency)
- Unique selectivity
- Straight-forward column packing

Go to bio-rad.com/CHTXT and explore the various features of CHT XT.

Ordering Information

Catalog #	Description
CHT Ceramic Hydroxyapatite XT Media	
12002457	CHT Ceramic Hydroxyapatite XT, 40 µm, 10 g
12002454	CHT Ceramic Hydroxyapatite XT, 40 µm, 100 g
12002456	CHT Ceramic Hydroxyapatite XT, 40 µm, 1 kg
12002455	CHT Ceramic Hydroxyapatite XT, 40 µm, 5 kg
Foresight Prepacked Columns	
12003150	Foresight CHT XT Column, 40 µm, 1 ml
12003149	Foresight CHT XT Column, 40 µm, 5 ml
Foresight Prepacked Chromatography Filter Plates*	
12003151	Foresight CHT XT Plates, 40 µm, 20 µl
Foresight RoboColumn Units**	
12003152	Foresight CHT XT RoboColumn Units, 40 µm, 200 µl
12003148	Foresight CHT XT RoboColumn Units, 40 µm, 600 µl

Services and Support

We offer multiple service and support levels to deliver solutions tailored to your needs.

- Method development
- Process development
- Custom resins
- Column packing support
- Global technical support



Process purification bottles, columns, and plates.

Global Supplier

We are a global supplier of process resins with more than 30 support sites across six continents. Contact your regional Bio-Rad Process Chromatography Specialist at process@bio-rad.com or call our customer service at 1-800-4-BIORAD (1-800-424-6723). Connect with us for the latest product information.



[linkedin.com/company/bio-rad](https://www.linkedin.com/company/bio-rad)



twitter.com/bioradlifesci

Screen Our Process Resins Today



Go to bio-rad.com/ResinSample and request a sample today.

Bio-Rad and CHT are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.

* Package size: 2 x 96-well plates.

** Package size: one row of eight columns.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 01 877 89019 **Belgium** 32 03 710 53 00 **Brazil** 55 11 3065 7550
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 36 01 459 6192 **Denmark** 45 04 452 10 00 **Finland** 35 08 980 422 00 **France** 33 01 479 593 00
Germany 49 089 3188 4393 **Hong Kong** 852 2789 3300 **Hungary** 36 01 459 6190 **India** 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 49486600
Japan 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 0 318 540 666 **New Zealand** 64 9 415 2280 **Norway** 47 0 233 841 30
Poland 36 01 459 6191 **Portugal** 351 21 4727717 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 36 01 459 6193 **Spain** 34 091 49 06 580
Sweden 46 08 555 127 00 **Switzerland** 41 0617 17 9555 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300
United Kingdom 44 01923 47 1301

