

# Got Purification Troubles?

## Minimize On-Column Aggregates during mAb Purification



### The Challenge

Have you seen two-peak elution of target monoclonal antibodies (mAb) after cation exchange (CEX) chromatography? The second elution peak consists of a mixture of monomers and aggregates of your target mAb and it leads to a decrease in your process yield.



### Purification Tip

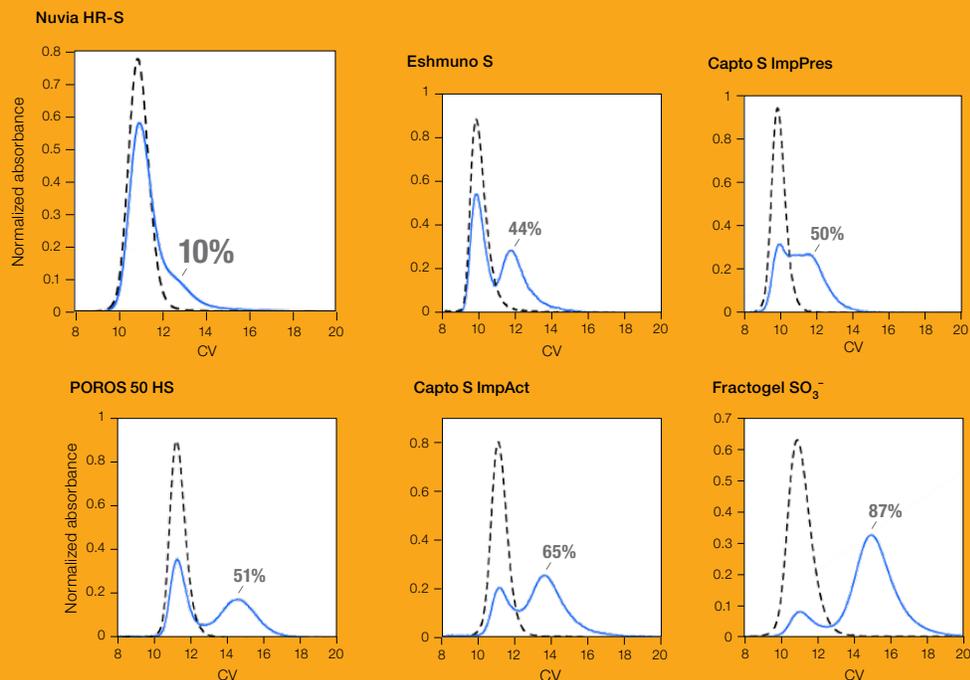
Choosing the optimal resin for your process will minimize on-column aggregates and maximize recovery. Be sure to screen multiple resins during your mAb purification process development.



### Our Approach

On-column aggregate formation depends on the specific chromatography resin used. Resin structure, surface extenders, pore size and distribution each contribute to on-column aggregate formation. Bio-Rad's Nuvia™ HR-S Resin — a strong cation exchanger with a pore size of 50  $\mu\text{m}$  without surface extenders — minimizes on-column aggregates. Nuvia HR-S generates only 10% aggregate formation relative to competing resins under similar experimental conditions.

**Fig. 1.** 0 min hold time (---); 1,000 min hold time (—); % of on-column aggregate shown in plots. For further details, refer to Bio-Rad bulletin 7000.



Go to [bio-rad.com/Nuvia-HR-S](http://bio-rad.com/Nuvia-HR-S) and explore Nuvia HR-S.



**BIO-RAD**

# Innovative Products. Customized Solutions.

## Nuvia™ HR-S Resin

Nuvia HR-S will put you on the path to:

- Sharper elution peaks
- Maximized process recovery
- Higher resolution between charge variants
- More robust process development

Go to [bio-rad.com/Nuvia-HR-S](http://bio-rad.com/Nuvia-HR-S) and explore Nuvia HR-S.

### Ordering Information

Catalog #	Description
732-4707	Foresight Nuvia HR-S Plates, 2 x 96-well, 20 µl
732-4831	Foresight Nuvia HR-S RoboColumn Unit, 200 µl
732-4832	Foresight Nuvia HR-S RoboColumn Unit, 600 µl
732-4723	Foresight Nuvia HR-S Column, 1 x 1 ml
732-4743	Foresight Nuvia HR-S Column, 1 x 5 ml
1560511	Nuvia HR-S Resin, 25 ml
1560513	Nuvia HR-S Resin, 100 ml
156-0515	Nuvia HR-S Resin, 500 ml
156-0517	Nuvia HR-S Resin, 10 L

## 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](mailto: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](https://twitter.com/bioradlifesci)

## Screen Our Process Resins Today



Go to [bio-rad.com/ResinSample](http://bio-rad.com/ResinSample) and request a sample today.

Capto is a trademark of General Electric Company. Fractogel and Eshmuno are trademarks of Merck KGaA, Darmstadt, Germany. POROS is a trademark of Thermo Fisher Scientific.

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [bio-rad.com](http://bio-rad.com) **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 1 877 89 01 177 **Belgium** 32 (0)3 710 53 00 **Brazil** 55 11 3065 7550 **Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 45 44 52 10 00 **Finland** 358 09 804 22 00 **France** 33 01 47 95 69 65 **Germany** 49 89 31 884 0 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 216091 **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 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 (0) 861 246 723 **Spain** 34 91 590 5200 **Sweden** 46 08 555 12700 **Switzerland** 41 026 674 55 05 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300 **United Kingdom** 44 020 8328 2000

