

Got Purification Troubles?

Get the Best HCP and Aggregate Clearance during mAb Purification



The Challenge

Are you unable to efficiently reduce process- and product-related impurities during your purification process? If you are performing flow-through purification, host cell protein (HCP) and aggregate removal could be poor in the flow-through fraction due to nonoptimal conditions, a nonideal resin, or both.



Purification Tip

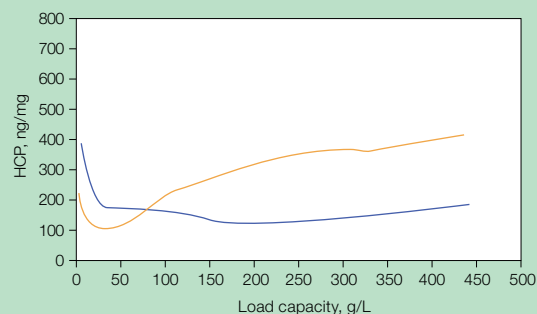
Screening multiple resins helps you select the right one for your purification success.



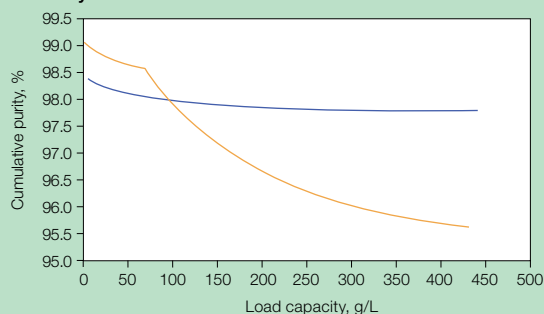
Our Approach

We tested the efficiency of two commercial hydrophobic anion exchange mixed-mode chromatography resins for the purification of an antibody intermediate (pI 9.12) from a Protein A resin eluate containing HCP levels and aggregate as indicated in Figures 1A and B. Nuvia aPrime 4A Resin consistently demonstrated the best reduction of HCP and aggregates.

A. HCP clearance



B. Purity



Resin	Remaining HCP Content, ng/mg	Purity, %
Nuvia aPrime 4A	159	97.8
Capto adhere	384	95.7

Fig. 1. Plots from the respective 1 ml columns with remaining HCP content indicated in the table at a loading concentration of 400 mg/ml. A, starting HCP content: 1,116 ng/mg. Flow-through conditions: Tris-based buffer system at pH 7.4. HCP content was measured by enzyme-linked immunosorbent assay. **B,** starting monomer: 94.8%. Purity was determined by ultra performance liquid chromatography. Figures 1A and B: Capto adhere Resin (—); Nuvia aPrime 4A Resin (—).

Data courtesy of AstraZeneca.

Go to [bio-rad.com/NuviaaPrime4A](https://www.bio-rad.com/NuviaaPrime4A) and explore the various features of Nuvia aPrime 4A Resin!



BIO-RAD

Innovative Products. Customized Solutions.

Nuvia aPrime 4A Resin

Nuvia aPrime 4A Resin delivers:

- A wide experimental design space
- Optimal product recovery
- Single-step removal of multiple impurities

Go to bio-rad.com/NuviaaPrime4A and explore the various features of Nuvia aPrime 4A Resin!

Ordering Information

Catalog #	Description
12007397	Nuvia aPrime 4A Resin, 25 ml
12007396	Nuvia aPrime 4A Resin, 100 ml
12007379	Nuvia aPrime 4A Resin, 500 ml
12007380	Nuvia aPrime 4A Resin, 5 L
12007391	Nuvia aPrime 4A Resin, 10 L
12007411	Foresight Nuvia aPrime 4A Plates, 20 µl
12007392	Foresight Nuvia aPrime 4A Column, 1 ml
12007393	Foresight Nuvia aPrime 4A Column, 5 ml
12007394	Foresight Nuvia aPrime 4A RoboColumn Unit, 200 µl
12007395	Foresight Nuvia aPrime 4A RoboColumn Unit, 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 is a trademark of Bio-Rad Laboratories, Inc. All trademarks used herein are the property of their respective owner.

Nuvia aPrime 4A Resin is covered by U.S. Patent Number 9,669,402 and foreign counterparts.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website bio-rad.com USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 00 800 00 24 67 23 Belgium 00 800 00 24 67 23 Brazil 4003 0399
Canada 1 905 364 3435 China 86 21 6169 8500 Czech Republic 00 800 00 24 67 23 Denmark 00 800 00 24 67 23 Finland 00 800 00 24 67 23
France 00 800 00 24 67 23 Germany 00 800 00 24 67 23 Hong Kong 852 2789 3300 Hungary 00 800 00 24 67 23 India 91 124 4029300 Israel 0 3 9636050
Italy 00 800 00 24 67 23 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Luxembourg 00 800 00 24 67 23 Mexico 52 555 488 7670
The Netherlands 00 800 00 24 67 23 New Zealand 64 9 415 2280 Norway 00 800 00 24 67 23 Poland 00 800 00 24 67 23 Portugal 00 800 00 24 67 23
Russian Federation 00 800 00 24 67 23 Singapore 65 6415 3188 South Africa 00 800 00 24 67 23 Spain 00 800 00 24 67 23 Sweden 00 800 00 24 67 23
Switzerland 00 800 00 24 67 23 Taiwan 886 2 2578 7189 Thailand 66 2 651 8311 United Arab Emirates 36 1 459 6150 United Kingdom 00 800 00 24 67 23

