

Got Purification Troubles?

Get the Best Monomer Recovery during Protein Purification



The Challenge

Is your protein unable to bind efficiently to your resin? If you are performing flow-through (FT) purification, monomer recovery could be poor in the flow-through fraction due to nonoptimal conditions, a nonideal resin, or both.



Purification Tip

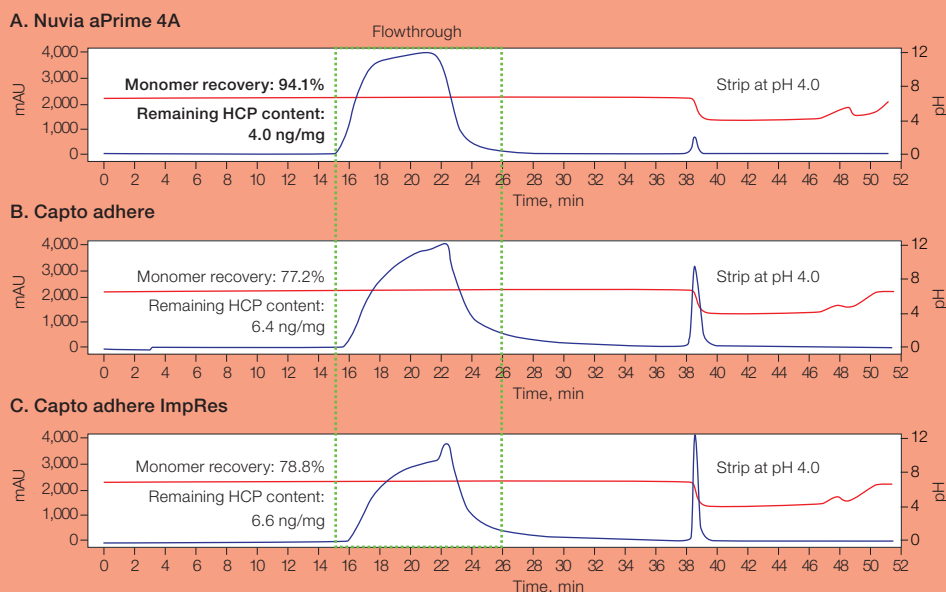
Performing a design of experiment (DOE) study of different pH, salt concentrations, and resins will help you select the best conditions and resin for your purification success.



Our Approach

A DOE study showed that Nuvia aPrime 4A Resin offers better monomer recovery than both Capto adhere Media for the purification of a basic protein (approx. pI of 8.4) at the optimal FT operation pH of 6.5. Tailing of the FT peak was observed from both Capto adhere Columns but not the Nuvia aPrime 4A Column. A spike at 280 nm was also seen when the Capto adhere Columns were washed with additional equilibration buffer for recovering target protein monomer. Nuvia aPrime 4A Resin's superior performance is the result of its optimized pore structures and surface properties, which offer efficient mass transfer and minimal nonspecific interaction between biomolecules and this chromatography media (Figure 1).

Fig. 1.
Chromatograms from the respective 1 ml columns with monomer recovery and remaining HCP content indicated. FT conditions: 99% monomer content in sample load, 150–200 ng/mg starting HCP content, 20 mM NaPi, 100 mM NaCl, pH 6.5. OD 280 (—); pH (—).



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



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!

Ordering Information

Catalog #	Description
12007397	Nuvia aPrime 4A, 25 ml
12007396	Nuvia aPrime 4A, 100 ml
12007379	Nuvia aPrime 4A, 500 ml
12007380	Nuvia aPrime 4A, 5 L
12007391	Nuvia aPrime 4A, 10 L
12007411	Foresight Nuvia aPrime 4A Plate, 20 μ l
12007392	Foresight Nuvia aPrime 4A Column, 1 ml
12007393	Foresight Nuvia aPrime 4A Column, 5 ml
12007394	Foresight Nuvia aPrime 4A RoboColumn, 200 μ l
12007395	Foresight Nuvia aPrime 4A RoboColumn, 600 μ 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 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. in certain jurisdictions. All trademarks used herein are the property of their respective owner.

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 4727117 **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

