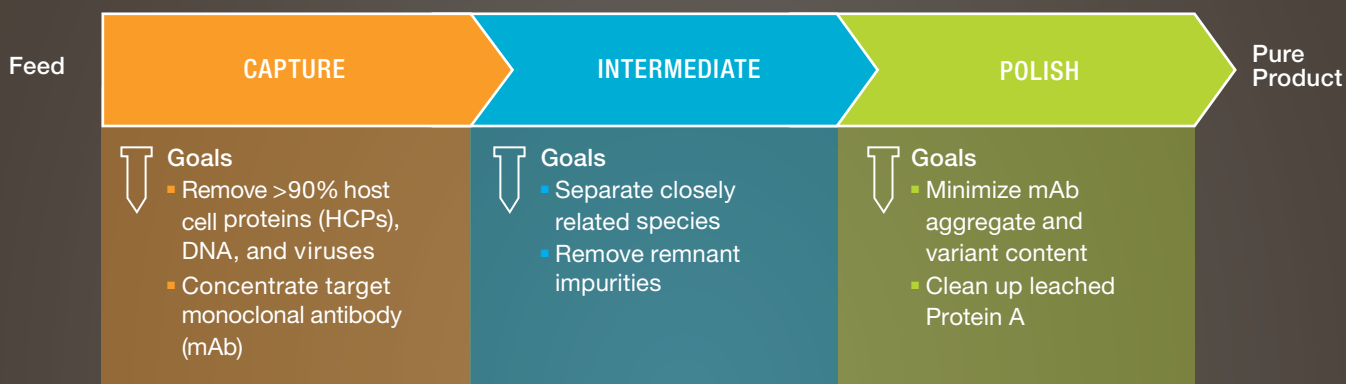
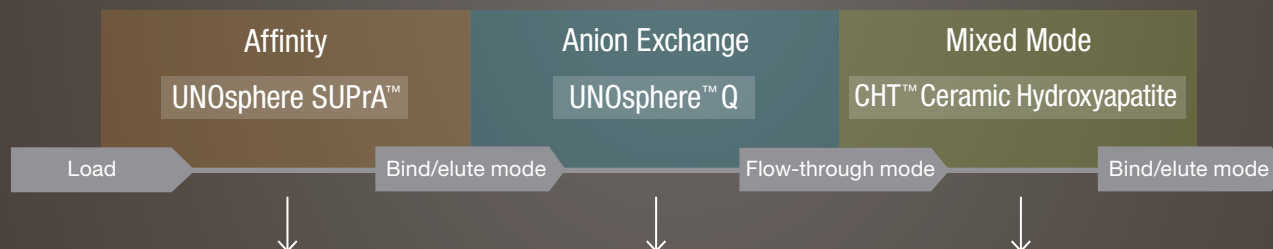




Quick Purification Strategy Affinity-Based Capture



Recommended Resins



Advantages

- Same set of buffers works for multiple mAbs (reduced optimization)
- Target molecule is in flowthrough, therefore less optimization needed relative to using cation exchange in bind/elute mode
- Best aggregate, endotoxin, and Protein A removal

Results

| | Feed | Capture | Intermediate | Polish |
|------------------|------------------------|---------|--------------|--------|
| Protein A, ng/mg | NA | ND | 112 | <0.4 |
| HCP, ng/mg | 1.4 x 10 ⁶ | 197 | 86 | 48 |
| DNA, ng/mg | >1.6 x 10 ⁵ | 19 | 1.9 | 3 |
| Aggregate % | ND | 42 | 40 | <0.1 |

NA, not applicable. ND, not determined.

Refer to experimental details on page 2.

Monoclonal Antibody Purification Strategy Details

Capture Conditions

| Step | Buffer | Column Volume | Flow Rate, cm/hr |
|----------------|------------------------|---------------|------------------|
| Equilibration | 1x PBS | 10 | 300 |
| Sample Loading | CHO cell supernatant | - | 300 |
| Wash | 1x PBS | 20 | 300 |
| Elution | 100 mM glycine, pH 3.0 | - | 300 |

Pooled elution fractions incubated for 1 hr at pH 3.0 to mimic virus deactivation.

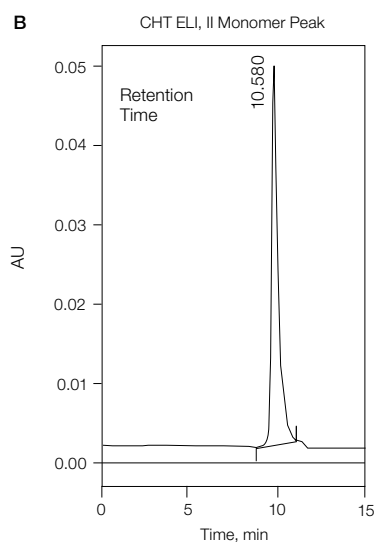
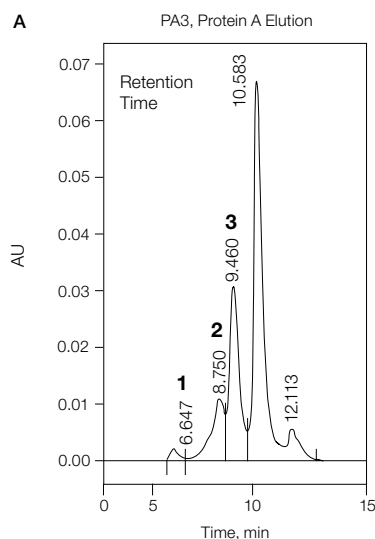
Intermediate Conditions

| Step | Buffer | Column Volume | Flow Rate, cm/hr |
|-------------------|---|---------------|------------------|
| Equilibration | 10 mM Na ₃ PO ₄ , pH 7.8 (buffer A) | 10 | 300 |
| Sample Loading | Eluate from UNOsphere SUPra column adjusted to pH 7.8 | - | 300 |
| Flow-Through Wash | Buffer A | 10 | 300 |

Flow-through fractions collected.

Polish Conditions

| Step | Buffer | Column Volume | Flow Rate, cm/hr |
|----------------|---|---------------|------------------|
| Equilibration | 10 mM Na ₃ PO ₄ , pH 6.8 (buffer A) | 10 | 300 |
| Sample Loading | UNOsphere Q flow-through pooled fractions adjusted to pH 6.8 | - | 300 |
| Wash | Buffer A | 15 | 300 |
| Elution | Salt gradient between buffer A and 10 mM Na ₃ PO ₄ , 1 M NaCl, pH 6.8 | 25 | 300 |



Removal of impurities by CHT chromatography. Elution fractions from the capture (**A**) and final polishing (**B**) chromatographic steps were analyzed by HPLC-SEC using a Bio-Sil® SEC 400-5 Column to evaluate levels of contamination. The peaks corresponding to aggregates (peaks 1 and 2) and dimers (peak 3) are removed by CHT to yield a clean sample. AU, absorbance units.

Resources

Download [bulletin 5728](#) to get details about this workflow.

Visit bio-rad.com/web/ProcessResins to get technical details about the resins used in this workflow.

Visit bio-rad.com/web/ProcessApplications to see other applications in which these resins can be used.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site 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 026674 55 05 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300 **United Kingdom** 44 020 8328 2000

