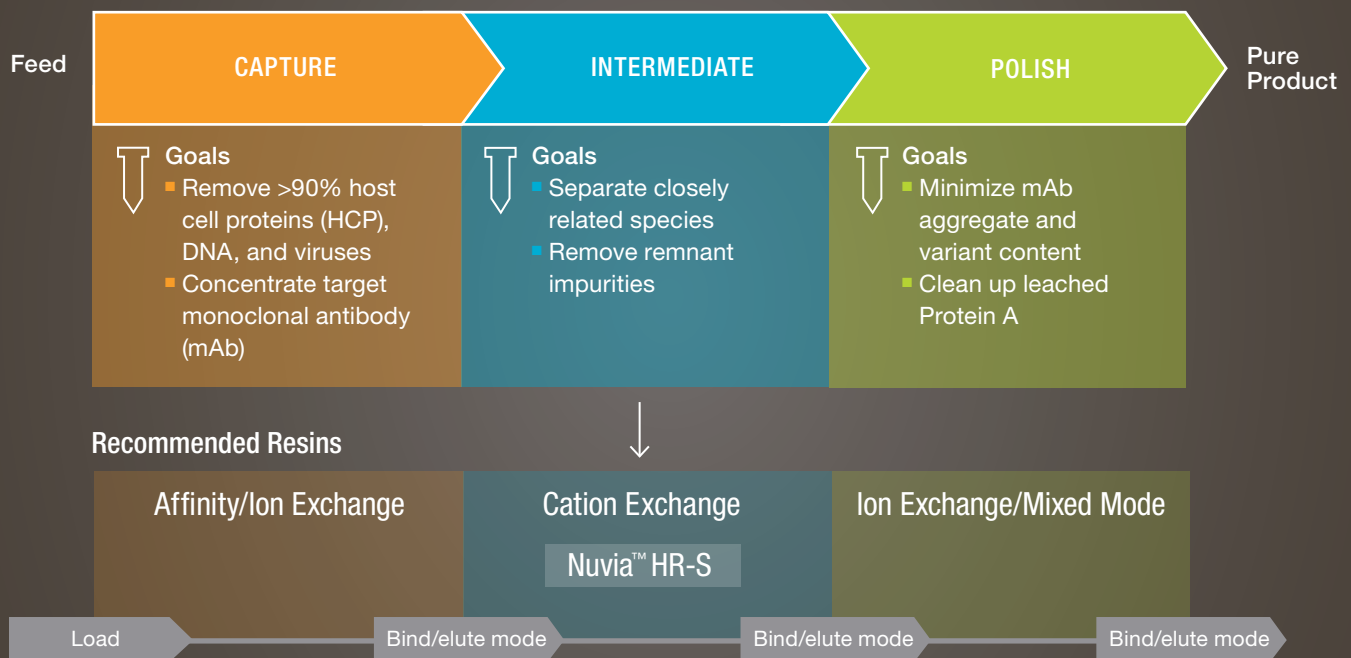




Quick Purification Strategy Resolving Monoclonal Antibody Charge Variants



Protein deamidation at neutral or alkaline pH is of significant concern to manufacturers of biotherapeutic proteins.

Deamidation at asparagine and glutamine residues:

Caused by

- Enzymatic or alkaline hydrolysis

Affects

- Immunogenicity, stability, and activity of proteins

Results in

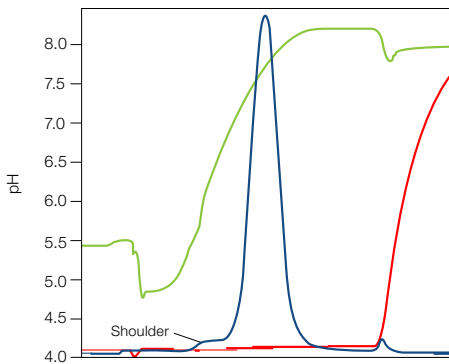
- Minimal changes in protein structure and charge
- Downstream purification challenges due to deaminated species
- Molecular heterogeneity

Refer to experimental details on page 2.

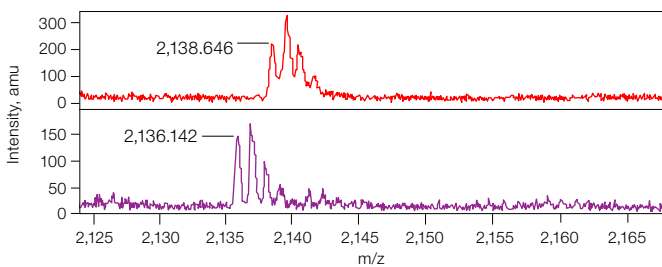
Monoclonal Antibody Purification Strategy Details

Intermediate purification to resolve charge variants with Nuvia HR-S

	Step	Buffer	Column Volume	Flow Rate, cm/hr	
Capture Affinity/ion exchange chromatography	Equilibration	10 mM sodium acetate, pH 4.5 (buffer A)	15	300	Polish Ion exchange/mixed-mode chromatography
	Sample Loading	Eluate from the capture step diluted 1:4 in buffer A	-	300	
	Wash 1	Buffer A	15	300	
	Wash 2	10 mM sodium citrate, 10 mM sodium phosphate, pH 5.0 (buffer B)	15	300	
	Elution	Gradient of buffer B to 10 mM sodium citrate, 10 mM sodium phosphate, pH 8.0	15	300	



Purification of the mAb on Nuvia HR-S. The distinct shoulder on the front side of the main peak is a modified mAb fragment. A₂₈₀ (—); conductivity (—); pH (—).



Mass spectrometry (MS) analysis of the mAb. MS analyses of the shouldered peak (top spectrum) and the main peak (lower spectrum) show a mass shift of about 2.5 atomic mass units between the two peaks, which is consistent with two or three deamidation events on a single peptide.

10	20	30	40	50	60	70
MYLLPATAA	GLLLAAQPA	HADIQTQSP	SSLASVGDV	VTITCRASQS	ISSYLNWYQQ	KPGKAPKLLI
80	90	100	110	120	130	140
YAASLQSGV	PSRFSQSGG	TDFTLTSSL	QPEDFATYYC	QQSYSTLLTF	GGGKVEIKR	TVAAPSVFIF
150	160	170	180	190	200	210
PPSDEQLRSG	TASVVCLLNN	FYPREVKVQV	K ¹⁸⁰ VDNALQSGNS	Q ¹⁸⁵ QESVTEQDSK	DSTYLSLST	LTLKADYEK
220	230	240	250	260		
HKVYACEVTH	QGLSSPVTKS	FNRGEYSARQ	STPPVCEYQG	QSSDLP		

VDNALQSGNSQESVTEQDSK

Shoulder fraction protein sequence. Examination of the shoulder fraction protein sequence confirms deamidation events on a single mAb peptide.

Resources

Visit bio-rad.com/web/NuviaHRS to get more information.

Visit bio-rad.com/web/ProcessResins to get technical details about Nuvia HR-S Resin.

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

