



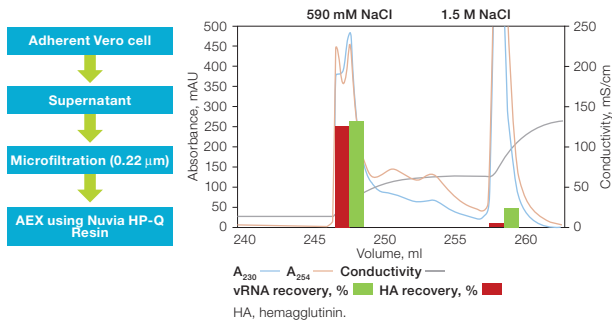
Quick Purification Strategy H1N1 Purification

Efficient virus recovery and impurity removal are challenges in the purification of viruses such as influenza A virus subtype H1N1. As a capture chromatography medium, Nuvia HP-Q Resin resulted in high virus recovery and excellent impurity removal in a single purification step. Moreover, Nuvia HP-Q Resin does not require benzonase treatment or the intensive filtration step of H1N1 culture harvest before loading. Here, we summarize H1N1 purification on Nuvia HP-Q Resin.

Procedure

H1N1 Purification on Nuvia HP-Q Resin

Format: Foresight Nuvia HP-Q 1 ml Column, 5.83 x 10¹¹ viral RNA (vRNA) copies/ml resin



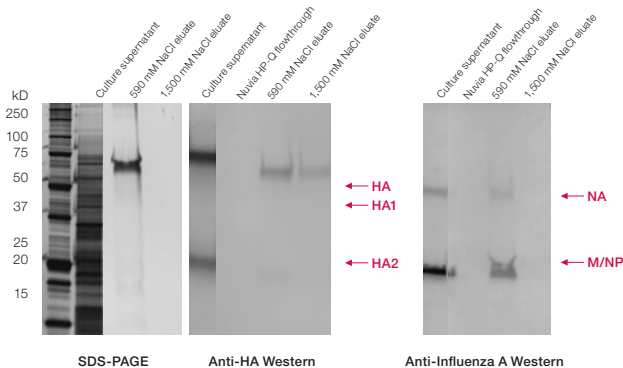
Optimized Protocol for Nuvia HP-Q Resin

Step	Mobile Phase	pH	CV	Remark
Equilibration	20 mM Tris, 100 mM NaCl	7.5	5	
Sample loading	20 mM Tris, 100 mM NaCl	7.5	-	5.83 x 10 ¹¹ viral RNA copies/ml of resin
Wash	20 mM Tris, 100 mM NaCl	7.5	10-30	
Elution	20 mM Tris, 590 mM NaCl	7.5	10	
High salt strip	20 mM Tris, 1.5 M NaCl	7.5	5	
Clean in place	1 M NaOH	>12	10	

* Operational linear flow rate: 84 cm/hr.

Results

SDS-PAGE and Western Blot Analyses



Average size: 128 ± 4.18 nm
Reported size range: 80-180 nm

Purified H1N1 virus was highly monodispersed with 99.8% monomer with a diameter of 128 nm.

Data generated by Bioprocessing Technology Institute (BTI) Singapore.

Results

Purification Step	HA Recovery	HCP/HA, mg/mg	Host cell DNA/HA, ng/mg
Culture supernatant	100.0	13.9	442.5
Microfiltration	100.0	13.0	221.3
Nuvia HP-Q eluate	80.0	1.1	0.3
EP guideline (for 15 HA dose)*	–	≤6.7	≤0.67

* ema.europa.eu/en/documents/scientific-guideline/guideline-influenza-vaccines-quality-module-revision-1_en.pdf

HA, hemagglutinin; HCP, host cell protein.

Highlights

- Benzonase treatment not required in this workflow
- High binding capacity
- Effective impurity removal

Reported Purification Schemes	HA Recovery, %	HCP/HA, mg/mg	hcDNA/HA, ng/mg	Reference, DOI
Diafiltration/microfiltration/ultrafiltration size exclusion chromatography, anion exchange chromatography (AEC)	53.0	7.7	34.5	10.1002/bit.21109
AEC, hydrophobic interaction chromatography	92.2	17.2	2.0	10.1016/j.jchromb.2019.03.037
AEC monolith	89.0	NA	NA	10.1016/j.jchromb.2017.06.086
Pseudo-affinity (saturated membrane)	94.2	2.1	23.0	10.1002/bit.22345
Pseudo-affinity (saturated membrane)	64.0	13	3.8	10.1002/jctb.5474

Visit bio-rad.com/NuviaHP-Q to get technical details about the resin used in this workflow.

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

Visit bio-rad.com/processlibrary to speak with a scientist or request a customized purification seminar.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc. All trademarks used herein are the property of their respective owner.



**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

