



Virus Purification Resin Selection Card

Bio-Rad offers a wide range of resins for each phase of virus process purification, at scales ranging from nanograms to kilograms.

Use this card to select the optimal resin for your virus process purification.



Visit bio-rad.com/web/VirusChromResins for more information.

Optimal Resin for Virus Purification

Property	Process Purification Resin Type					
	Adenovirus			Other Small to Midsized Mammalian Viruses		
Resin type	Nuvia™ cPrime™	Nuvia Q	UNOsphere™ Q	CHT™ Ceramic Hydroxyapatite	CFT™ Ceramic Fluoroapatite	MPC™ Ceramic Hydroxyfluoroapatite
Chromatography type	Mixed-mode (HIC and cation exchange)	Anion exchange	Anion exchange	Mixed-mode (cation exchange and metal affinity)	Mixed-mode (cation exchange and metal affinity)	Mixed-mode (cation exchange and metal affinity)
Particle size	70 ± 10 µm	85 ± 15 µm	120 µm	20 ± 2, 40 ± 4, 80 ± 8 µm	40 ± 4 µm	40 ± 4 µm
Ionic capacity	110–150 µeq/ml	100–170 µeq/ml	120 µeq/ml	—	—	—
Dynamic binding capacity (DBC)	>40 mg hlgG/ml (at 10% breakthrough) at 300 cm/hr* >60 mg lactoferrin/ml*	≥170 mg/ml at 300 cm/hr*	≥180 mg BSA/ml at 150 cm/hr* ≥125 mg BSA/ml at 600 cm/hr*	≥25 mg lysozyme/g* 25–60 mg IgG/ml at 300 cm/hr*	14–21 mg lysozyme/g* 33 mg IgG/ml*	≥25 mg lysozyme/g* 25–50 mg IgG/ml*
Recommended linear flow rate	50–600 cm/hr	50–600 cm/hr	50–300 cm/hr	50–300 cm/hr	50–300 cm/hr	50–300 cm/hr
pH stability	Short-term: 3–14 Long-term: 4–13	Short-term: 2–14 Long-term: 4–12	1–14	6.5–14	Operating pH: 5–14 Storage pH: 11–14	6.5–14

BSA, bovine serum albumin; HIC, hydrophobic interaction chromatography.

Refer to bulletins 6790 and 6807 to see how these resins help in virus process purification.

* Go to the product detail page or bulletin 6713 on bio-rad.com to see how the DBC was determined and for other technical details.

Ordering Information

Nuvia cPrime	Catalog #	Size	Nuvia Q	Catalog #	Size	UNOsphere Q	Catalog #	Size
Foresight™ Nuvia cPrime Plates	732-4705	2 x 96-well, 20 µl	Foresight Nuvia Q Plates	732-4703	2 x 96-well, 20 µl	Foresight UNOsphere Q Plates	732-4714	2 x 96-well, 20 µl
Foresight Nuvia cPrime RoboColumn Unit	732-4807 732-4808	200 µl 600 µl	Foresight Nuvia Q RoboColumn Unit	732-4804 732-4805	200 µl 600 µl	Foresight UNOsphere Q RoboColumn Unit	732-4819 732-4820	200 µl 600 µl
Foresight Nuvia cPrime Column	732-4722 732-4742	1 ml 5 ml	Foresight Nuvia Q Column	732-4721 732-4741	1 x 1 ml 1 x 5 ml	Foresight UNOsphere Q Column	732-4732 732-4752	1 ml 5 ml
Nuvia cPrime Resin	1563401 1563402 156-3403 156-3404 156-3405 156-3406	25 ml 100 ml 500 ml 1 L 5 L 10 L	Nuvia Q Resin	1560411 1560413 156-0415 156-0417	25 ml 100 ml 500 ml 10 L	UNOsphere Q Resin	1560101 1560103 156-0105 156-0107	25 ml 100 ml 500 ml 10 L

CHT Ceramic Hydroxyapatite	Catalog #	Size	CFT Ceramic Fluoroapatite	Catalog #	Size	MPC Ceramic Hydroxyfluoroapatite	Catalog #	Size
Foresight CHT Type I Plates	732-4716	40 µm, 2 x 96-well, 20 µl	Bio-Scale™ Mini CFT Type II Cartridges	7324405 7324406	1 x 5 ml 5 x 5 ml	Foresight MPC Type I Plates	732-4785	40 µm, 2 x 96-well, 20 µl
Foresight CHT Type II Plates	732-4718	40 µm, 2 x 96-well, 20 µl	CFT Ceramic Fluoroapatite Type II Media, 40 µm	1585200 1575000 157-5100 157-5500	10 g 100 g 1 kg 5 kg	Foresight MPC Type I RoboColumn Unit	732-4828 732-4829	40 µm, 200 µl 40 µm, 600 µl
Foresight CHT Type I RoboColumn Unit	732-4822	40 µm, 200 µl				Foresight MPC Type I Column	732-4737 732-4757	40 µm, 1 ml 40 µm, 5 ml
Foresight CHT Type II RoboColumn Unit	732-4823 732-4825 732-4826	40 µm, 600 µl 40 µm, 200 µl 40 µm, 600 µl				MPC Ceramic Hydroxyfluoroapatite Type I Media, 40 µm	1580200 1570200 157-0201 157-0205	10 g 100 g 1 kg 5 kg
Foresight CHT Type I Column	732-4735 732-4755	40 µm, 1 ml 40 µm, 5 ml						
Foresight CHT Type II Column	732-4736 732-4756	40 µm, 1 ml 40 µm, 5 ml						
CHT Type I Media, 40 µm	1584000 1570040 157-0041 157-0045	10 g 100 g 1 kg 5 kg						
CHT Type I Media, 80 µm	1588000 1570080 157-0081 157-0085	10 g 100 g 1 kg 5 kg						
CHT Type II Media, 40 µm	1584200 1574000 157-4100 157-4500	10 g 100 g 1 kg 5 kg						
CHT Type II Media, 80 µm	1588200 1578000 157-8100 157-8500	10 g 100 g 1 kg 5 kg						

RoboColumn is a trademark of Atoll GmbH.

