



## Aggregate Purification Resin Selection Card

Bio-Rad offers a wide range of resins for aggregate depletion during process-scale purification.

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



Visit [bio-rad.com/web/AggChromResins](http://bio-rad.com/web/AggChromResins) for more information.

### Optimal Resin for Aggregate Purification

Property	Process Purification Resin Type					
Resin type	Nuvia™ S	Nuvia HR-S	Macro-Prep® Methyl	Macro-Prep t-Butyl	CHT™ Ceramic Hydroxyapatite	Nuvia™ cPrime™
Chromatography type	Cation exchange	Cation exchange	Hydrophobic interaction	Hydrophobic interaction	Mixed-mode (cation exchange and metal affinity)	Mixed-mode (HIC and cation exchange)
Particle size	85 ± 15 µm	50 ± 10 µm	50 µm	50 µm	20 ± 2, 40 ± 4, 80 ± 8 µm	70 ± 10 µm
Ionic capacity	90–150 µeq/ml	100–180 µeq/ml	<2 µeq/ml	120 µeq/ml	—	110–150 µeq/ml
Dynamic binding capacity (DBC)	≥110 mg/ml at 300 cm/hr*	≥70 mg/ml at 300 cm/hr*	15 mg HSA/ml*	25 mg HSA/ml*	≥25 mg lysozyme/g* 25–60 mg IgG/ml at 300 cm/hr*	>40 mg hlgG/ml (at 10% breakthrough) at 300 cm/hr* >60 mg lactoferrin/ml*
Recommended linear flow rate	50–300 cm/hr	50–200 cm/hr	100–600 cm/hr	100–600 cm/hr	50–300 cm/hr	50–600 cm/hr
pH stability	Short-term: 2–14 Long-term: 4–13	Short-term: 2–14 Long-term: 4–13	1–10	1–10	6.5–14	Short-term: 3–14 Long-term: 4–13

HIC, hydrophobic interaction chromatography; HSA, human serum albumin.

Refer to [bulletin 6808](#) to see how these resins help in aggregate process purification.

\* Go to the product detail page or [bulletin 6713](#) on [bio-rad.com](http://bio-rad.com) to see how the DBC was determined and for other technical details.

### Ordering Information

Nuvia S	Catalog #	Size
Foresight™ Nuvia S Plates	732-4701	2 x 96-well, 20 µl
Foresight Nuvia S RoboColumn Unit	732-4801	200 µl
	732-4802	600 µl
Foresight Nuvia S Column	732-4720	1 x 1 ml
	732-4740	1 x 5 ml
Nuvia S Resin	1560311	25 ml
	1560313	100 ml
	156-0315	500 ml
	156-0317	10 L

Nuvia HR-S	Catalog #	Size
Foresight Nuvia HR-S Plates	732-4707	2 x 96-well, 20 µl
Foresight Nuvia HR-S RoboColumn Unit	732-4831	200 µl
	732-4832	600 µl
Foresight Nuvia HR-S Column	732-4723	1 x 1 ml
	732-4743	1 x 5 ml
Nuvia HR-S Resin	1560511	25 ml
	1560513	100 ml
	156-0515	500 ml
	156-0517	10 L

Macro-Prep Methyl	Catalog #	Size
Macro-Prep Methyl HIC Resin	1580080	25 ml
	1560080	100 ml
	156-0081	500 ml
	156-0082	5 L
	156-0083	10 L

Macro-Prep t-Butyl	Catalog #	Size
Macro-Prep t-Butyl HIC Resin	1580090	25 ml
	1560090	100 ml
	156-0091	500 ml
	156-0093	10 L

CHT Ceramic Hydroxyapatite	Catalog #	Size
Foresight CHT Type I Plates	732-4716	40 µm, 2 x 96-well, 20 µl
Foresight CHT Type II Plates	732-4718	40 µm, 2 x 96-well, 20 µl
Foresight CHT Type I RoboColumn Unit	732-4822	40 µm, 200 µl
	732-4823	40 µm, 600 µl
Foresight CHT Type II RoboColumn Unit	732-4825	40 µm, 200 µl
	732-4826	40 µm, 600 µl
Foresight CHT Type I Column	732-4735	40 µm, 1 ml
	732-4755	40 µm, 5 ml
Foresight CHT Type II Column	732-4736	40 µm, 1 ml
	732-4756	40 µm, 5 ml
CHT Type I Media, 40 µm	1584000	10 g
	1570040	100 g
	157-0041	1 kg
	157-0045	5 kg
CHT Type I Media, 80 µm	1588000	10 g
	1570080	100 g
	157-0081	1 kg
	157-0085	5 kg
CHT Type II Media, 40 µm	1584200	10 g
	1574000	100 g
	157-4100	1 kg
	157-4500	5 kg
CHT Type II Media, 80 µm	1588200	10 g
	1578000	100 g
	157-8100	1 kg
	157-8500	5 kg

Nuvia cPrime	Catalog #	Size
Foresight Nuvia cPrime Plates	732-4705	2 x 96-well, 20 µl
Foresight Nuvia cPrime RoboColumn Unit	732-4807	200 µl
	732-4808	600 µl
Foresight Nuvia cPrime Column	732-4722	1 ml
	732-4742	5 ml
Nuvia cPrime Resin	1563401	25 ml
	1563402	100 ml
	156-3403	500 ml
	156-3404	1 L
	156-3405	5 L
	156-3406	10 L

RoboColumn is a trademark of Atoll GmbH.

