

Bio-Rad: Your Trusted Partner for Process Chromatography Resins

Innovative Resins, Customized Solutions

Since 1957, Bio-Rad Laboratories, Inc. has been a leading global provider of life science and clinical diagnostic products. Bio-Rad provides chromatography resins for biopharmaceutical workflows, including resin screening, preparative and analytical uses, and biomanufacturing. Our resins are used in commercial settings by customers worldwide for their downstream bioprocessing. We continue to expand solutions to support each phase of the biotherapeutic purification process. Bio-Rad is compliant with FDA, EMA, CFDA, and PMDA guidelines and is ISO-certified.

Resins in Multiple Formats to Serve Your Process Purification Needs

Ion Exchange	Affinity	Mixed-Mode
Nuvia HR-S	Nuvia IMAC	Nuvia aPrime 4A
Nuvia S	Affi-Gel™ 10	Nuvia cPrime
Nuvia HP-Q	Affi-Gel Boronate	CHT™ Ceramic Hydroxyapatite I
Nuvia Q	Affi-Gel Blue	CHT Ceramic Hydroxyapatite II
Macro-Prep™ High S	CM Affi-Gel Blue	CHT Ceramic Hydroxyapatite XT
Macro-Prep High Q	DEAE Affi-Gel Blue	Bio-Gel™ HT
Macro-Prep High Q-3HT	Affi-Gel Protein A	Bio-Gel HTP
Macro-Prep CM	Affi-Prep™ Protein A	
Macro-Prep DEAE		
UNOsphere S		
UNOsphere Q		
Hydrophobic Interaction	Size Exclusion	Specialty
Macro-Prep Methyl	Bio-Gel A	AG™ Resins
Macro-Prep t-Butyl	Bio-Gel P	Chelex™ 100
Bio-Beads SM-2	Bio-Beads S-X	Aminex™ Resins
		Bio-Rex Resins



Resins are available in multiple formats including: Foresight™ Columns, Plates, and RoboColumn Units for screening; process-scale prepacked Foresight Pro Chromatography Columns, and in bulk.

Supporting Your Purification Success — from Development to Commercialization

Knowledge Tools



Library

Find complete product information, technical videos, application notes, and guides for all process chromatography resins.



Webinars

Register to attend upcoming webinars or view Ask the Expert series webcasts for best practices and strategies for performing successful purifications.



Events

Join us at an upcoming event. Meet our scientists and learn about purification solutions for your evolving research needs.

Resin Samples



Bio-Rad manufactures a wide range of chromatography resins for process separations, providing process developers with great flexibility and high productivity.



Request a customized on-site or virtual seminar with our process scientists.

Expert Advice



Our team of scientists is here to accelerate your method and process development for biopharmaceutical manufacturing.

Visit bio-rad.com/ProcessChrom for more information.

BIO-RAD, AFFI-GEL, AFFI-PREP, AG, AMINEX, BIO-GEL, CHELEX, CHT, FORESIGHT, and MACRO-PREP are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner. © 2024 Bio-Rad Laboratories, Inc.

Nuvia aPrime 4A Resin is covered by U.S. Patent Number 9,669,402 and foreign counterparts.



**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** 82 080 007 7373 **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

