



Bio-Rad: Your Trusted Partner for Process Chromatography Resins

Innovative Products, Customized Solutions

Bio-Rad, a leading global provider of life science and clinical diagnostic products for more than 60 years, is an established supplier of process chromatography resins that provides researchers with the tools they need for isolating and purifying proteins from laboratory scale to bioprocess manufacturing. Our resins are used in many commercial settings and we continue to expand solutions to support each phase of the biotherapeutic purification process. Bio-Rad is FDA, EMA, CFDA, and PMDA compliant and ISO certified.

Resins in Multiple Formats to Serve Your Process Purification Needs

Ion Exchange

Nuvia HR-S
 Nuvia Q
 Nuvia S
 Nuvia HP-Q
 Macro-Prep CM
 Macro-Prep DEAE
 Macro-Prep High Q
 Macro-Prep High Q-3HT
 Macro-Prep High S
 UNOsphere Q
 UNOsphere S

Affinity

Nuvia IMAC
 UNOsphere SUPrA
 Affi-Gel Blue
 CM Affi-Gel Blue
 DEAE Affi-Gel Blue
 Affi-Gel Protein A
 Affi-Prep Protein A

Mixed Mode

Nuvia cPrime
 Nuvia aPrime 4A
 CHT Ceramic Hydroxyapatite I/II
 CHT Ceramic Hydroxyapatite XT
 MPC Ceramic Hydroxyfluoroapatite
 CFT Ceramic Fluoroapatite

Hydrophobic Interaction

Macro-Prep Methyl
 Macro-Prep t-Butyl

Size Exclusion

Bio-Gel P
 Bio-Gel A
 Bio-Beads SM-2
 Bio-Beads S-X

Specialty

AG Resins
 Chelex 100



Bottles and Foresight Columns, Plates, and RoboColumn Units

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.

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

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

BIO-RAD

**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

