



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
 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™
 CHT™ Ceramic Hydroxyapatite
 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

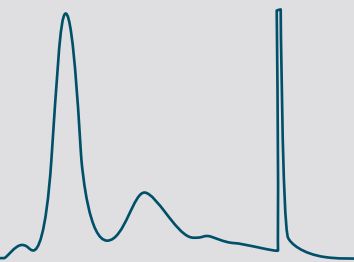
AG® Resins
 Chelex® 100



Bottles and Foresight™ Columns, Plates, and RoboColumn Units

Process Resin Services

- Methods development
- Process development
- Custom process resins
- Column packing support
- Global technical support



Process Purification Workflow

Bio-Rad's resins are used for multiple applications at different steps in the process purification workflow.

Capture

Intermediate

Polish

- Monoclonal antibodies
- Therapeutic enzymes, proteins, and hormones
- Viruses and vaccines
- Blood products
- Radio embolization
- Aggregate removal
- Endotoxin purification
- Biosimilars/biobetters
- And more...

Global Supplier

We are a global supplier of process resins with more than 30 support sites across six continents. Contact your regional Bio-Rad Process Chromatography Specialist at process@bio-rad.com or call our customer service at 1-800-4-BIORAD (1-800-424-6723). Connect with us for the latest product information.



[linkedin.com/company/bio-rad](https://www.linkedin.com/company/bio-rad)



[facebook.com/biorad](https://www.facebook.com/biorad)



twitter.com/bioradlifesci



[youtube.com/user/BioRadLifeScience](https://www.youtube.com/user/BioRadLifeScience)

Screen Our Process Resins Today

Go to [bio-rad.com/ResinSample](https://www.bio-rad.com/ResinSample) to request a sample.



RoboColumn is a trademark of Atoll GmbH.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 1 877 89 01 177 **Belgium** 32 (0)3 710 53 00 **Brazil** 55 11 3065 7550
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 45 44 52 10 00 **Finland** 358 09 804 22 00
France 33 01 47 95 69 65 **Germany** 49 89 31 884 0 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 972 03 963 6050 **Italy** 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 (0)318 540 666
New Zealand 64 9 415 2280 **Norway** 47 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04
Singapore 65 6415 3188 **South Africa** 27 (0) 861 246 723 **Spain** 34 91 590 5200 **Sweden** 46 08 555 12700 **Switzerland** 41 026 674 55 05
Taiwan 886 2 2578 7189 **Thailand** 66 662 651 8311 **United Arab Emirates** 971 4 8187300 **United Kingdom** 44 020 8328 2000

