



# 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

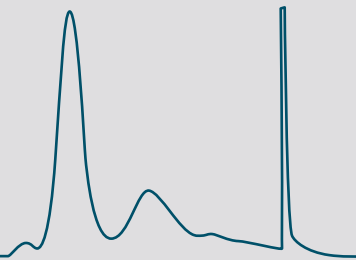
AG® Resins  
 Chelex® 100



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](mailto: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](https://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](http://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

