



## Bio-Rex™ Resins

### Family Members

- Bio-Rex 5 Resin
- Bio-Rex 70 Resin
- Bio-Rex Reactor Grade Resin

### At a Glance

The Bio-Rex Resins are weak ion exchangers. They are available in various sizes to aid laboratory- to manufacture-scale purifications.

### Bio-Rex 5 Resins

#### Features

- Weak anion exchangers that contain mostly tertiary and some quaternary amines
- Exhaustively sized, purified, and converted to make them suitable for accurate, reproducible analytical techniques
- Available in two dry mesh sizes (100–200 and 200–400) with analytical grade purity

#### Bio-Rex 5 Technical Specifications

Functional group	R-N <sup>+</sup> (CH <sub>3</sub> ) <sub>2</sub> C <sub>2</sub> H <sub>4</sub> OH
Matrix	Styrene divinylbenzene
Crosslinkage	4%
Approx. MW exclusion	1,400
Operating stability/pH range	<7/0–14
Min. wet capacity (meq/ml)	1.1
Nominal density	0.7
Ionic form	Chloride

#### Applications

- Have an affinity for organic acids that can be used to remove acids from sugars
- Exchange anions such as iodide, bromide, and chloride with weak interactions

### Bio-Rex 70 Resins

#### Features

- Weak cation exchangers that contain mostly tertiary and some quaternary amines
- Functional group: carboxylic acid exchange groups on a macroreticular acrylic polymer lattice
- Available in both analytical and biotechnology grades
  - Analytical grade: exhaustively sized, purified, and converted to make it suitable for accurate, reproducible analytical techniques
  - Biotechnology grade: analytical grade resin that has been certified to contain less than 100 microorganisms per gram of resin

#### Applications

- Purification and fractionation of peptides, proteins, antibiotics, and other cationic molecules

## Bio-Rex Reactor Grade Technical Specifications

Resin	Crosslinkage	Approx. MW Exclusion	Operating Stability/pH Range	Min. Wet Capacity	Nominal Density	Ionic Forms
<b>Bio-Rex MSZ 501(D) Resin</b>						
Type: Mixed bed	8 and 10%	1,000	0–14/0–14	1.5 meq/ml	0.72 g/ml	H <sup>+</sup> + OH <sup>-</sup>
Functional group: R-SO <sub>3</sub> <sup>-</sup> and R-CH <sub>2</sub> N <sup>+</sup> (CH <sub>3</sub> ) <sub>3</sub>						

## Bio-Rex Reactor Grade Resin

### Features

- Uniform bead size, which minimizes pressure drops and increases resin capacity; used for desalting water and other nonelectrolytes
- Contains a dye that changes from blue-green to gold when its exchange capacity is exhausted

### Applications

- Bio-Rex RG 501-X8 and Bio-Rex MSZ 501(D) Resins used for power plant deionization systems and large-scale cleanup of metals from wastewater

### Additional Information

Usage of Bio-Rex 5 Resin, [bulletin LIT207](#)

Usage of Bio-Rex 70 Resin, [bulletin LIT206](#)

Usage of Bio-Rex MSZ 501(D) Resin, [bulletin LIT205](#)

Large bulk volumes and special packaging for industrial applications are available on request.

For technical/product support or to request a quote, email your regional Bio-Rad representative or contact customer service at 1-800-4-BIORAD (1-800-424-6723).

Keep up with our updates by connecting with us.

 [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)

Visit [bio-rad.com/web/BioRexBeads](https://www.bio-rad.com/web/BioRexBeads) for more information and download our Process Resin Selection Guide ([bulletin 6713](#)) for technical and ordering information.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [bio-rad.com](https://www.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 2 651 8311 United Arab Emirates 971 4 8187300 United Kingdom 44 020 8328 2000

