



Bio-Gel® Resins

Family Members

- Bio-Gel P Polyacrylamide Gel
- Bio-Gel A Agarose Gel

At a Glance

Bio-Gel Size Exclusion Resins are engineered for gentle, high-resolution separation of sensitive compounds in aqueous or organic solutions. They are available in various mesh sizes to accommodate a wide range of applications. They can be stored at ambient temperatures and are available in various sizes to aid multiple scales of purification.

Bio-Gel P Technical Specifications

	Hydrated Bead Size, μm	Hydrated Bed Volume Mass Dry Gel, ml/g	Typical Flow Rate, cm/hr*	Molecular Weight Fractionation Range
Bio-Gel P-2, fine	45–90	3	5–10	100–1,800
Bio-Gel P-2, extra fine	<45		<10	
Bio-Gel P-4, medium	90–180	4	15–20	800–4,000
Bio-Gel P-4, fine	45–90		10–15	
Bio-Gel P-4, extra fine	<45		<10	
Bio-Gel P-6, medium	90–180	6.5	15–20	1,000–6,000
Bio-Gel P-6, fine	45–90		10–15	
Bio-Gel P-6, extra fine	<45		<10	
Bio-Gel P-6DG Gel	90–180		15–20	
Bio-Gel P-10 Gel, medium	90–180	7.5	15–20	1,500–20,000
Bio-Gel P-10 Gel, fine	45–90		10–15	
Bio-Gel P-30 Gel, medium	90–180	9	7–13	2,400–40,000
Bio-Gel P-30 Gel, fine	45–90		6–11	
Bio-Gel P-60 Gel, medium	90–180	11	4–6	3,000–60,000
Bio-Gel P-60 Gel, fine	45–90		3–5	
Bio-Gel P-100 Gel, medium	90–180	12	4–6	5,000–100,000
Bio-Gel P-100 Gel, fine	45–90		3–5	

* Flow rates determined in a 1.5 x 70 cm column using hydrostatic pressure head-to-bed ratio of 1:1

Bio-Gel P Polyacrylamide Gel

Features

- Prepared by copolymerization of acrylamide and *N,N'*-methylenebisacrylamide
- Chromatography type: size exclusion chromatography (SEC)
- Available in several particle size ranges with molecular weight exclusion limits ranging from 100 to 100,000 daltons
- Extremely hydrophilic and essentially nonionic
- Do not support microbial growth or leach carbohydrates (due to their synthetic composition), as dextrose and agarose gels can
- Autoclavable at pH 5.5–6.5 and operate over a pH range of 2–10
- Flow rate and resolution increase with increasing temperature in the range of 4–80°C

Applications

- Purification of nucleotides, peptides, and tannins

Related Product

Bio-Gel P-6DG Gel ([bulletin 2068](#)), with a molecular weight exclusion limit of 6,000 daltons, is recommended specifically for desalting and buffer exchange applications. It provides rapid results with high sample recovery and minimal sample dilution. Its hydrophilicity ensures little or no interaction between the gel and sample molecules.

Bio-Gel A Agarose Gel

Features

- Consists of agarose beads in which the pore size is controlled by the percentage of agarose in the gel
- Compatible with all commonly used buffers and can be used with high-salt buffers without significantly changing the bed volume
- Bio-Gel A 1.5 m gel may be used at pH 4–13 and at 2–30°C
- The fractionation range is from 10,000 to 1,500,000 daltons

Applications

- Purification of antibodies and aggregates
- Isolation of immune complexes
- DNA fractionation
- Chromatin isolation and purification

Additional Information

Usage of Bio-Gel P Polyacrylamide Gel, [bulletin LIT174](#)

Usage of Bio-Gel A Agarose Gel, [bulletin 4006139](#)

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

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

Visit [bio-rad.com/web/BioGelBeads](https://www.bio-rad.com/web/BioGelBeads) 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

