

Technical Bulletin

Gel Filtration Markers Kit for Protein Molecular Weights 29,000-700,000 Da

MWGF1000

Product Description

Gel filtration chromatography is an established method for determining the size and molecular mass of proteins. Fractionation is based on the diffusion of molecules into the pores of the resin. Larger proteins do not enter the pores of the resin as readily, but instead pass through the fluid volume of the column faster than smaller proteins. These protein molecules elute from the column in order of decreasing molecular mass.

The molecular mass determination of unknown proteins is made by comparing the ratio of V_e/V_o for the protein in question to the V_e/V_o of protein standards of known molecular mass, where:

- V_e is the elution volume
- V_o is the void volume

The V_o of a given column is based on the volume of effluent required for the elution of a large molecule such as Blue Dextran (molecular mass of ~2,000 kDa, Cat. No. D4772). Plotting the logarithms of the known molecular masses of protein standards versus their respective V_e/V_o values produces a linear calibration curve (such as shown in **Figure 1**).

V_e/V_o is essentially independent of column size and protein concentration. However, V_e/V_o may be dependent on temperature for some proteins. Unreliable molecular masses may be obtained if the protein:

- Forms a complex with the gel
- Contains a large amount of carbohydrate
- Aggregates to larger complexes
- Or dissociates into subunits under the conditions used.¹

The molecular mass of an impure protein may be determined using this procedure if a specific detection test is available for the protein.

The procedure for determining molecular masses using gel filtration chromatography, as outlined in this bulletin, is a modification of published methods.^{1,2} The protein standards in this kit may be suitable for use in other chromatographic systems such as HPLC, although some buffer systems seem to alter the elution volumes of albumin (Cat. No. A8531) and carbonic anhydrase (Cat. No. C7025). The proteins in the MWGF1000 Kit have a range of molecular masses from 29 kDa to 699 kDa.

Several theses³ and dissertations⁴⁻²⁷ cite use of product MWGF1000 in their protocols.

Reagent

- Carbonic Anhydrase from bovine erythrocytes, Cat. No. C7025, 15 mg/vial. Approximate Molecular Mass: ~29 kDa
- Albumin, Bovine Serum, Cat. No. A8531, 50 mg/vial (contains ~0.3% dithiothreitol). Approximate Molecular Mass: ~66 kDa
- Alcohol Dehydrogenase from yeast, Cat. No. A8656, 25 mg/vial. Approximate Molecular Mass: ~150 kDa
- β -Amylase from sweet potato, Cat. No. A8781, 15 mg protein/vial (contains ~15% NaCl, 4% glucose, and 1% dithiothreitol). Approximate Molecular Mass: ~200 kDa
- Apoferritin from horse spleen, Cat. No. A3660, 50 mg/vial (supplied as a solution at 25 mg protein/mL in 50% glycerol with 0.075 M NaCl). Approximate Molecular Mass: ~443 kDa
- Thyroglobulin, bovine, Cat. No. T9145, 35 mg/vial. Approximate Molecular Mass: ~669 kDa
- Blue Dextran, Cat. No. D4772, 50 mg/vial. Approximate Molecular Mass: ~2,000 kDa

Storage/Stability

Store the MWGF1000 product at -20 °C.

MILLIPORE
SIGMA

Procedure

Use of MWGF1000 for Gel Filtration Chromatography

Buffer and Resin:

- It is recommended to use 50 mM Tris-HCl, pH 7.5, with 100 mM KCl as the equilibration buffer, with a 90 cm × 1.6 cm Sepharose® CL-6B (Cat. No. CL6B200) column at 2-8 °C.
- For information on resin preparation, column packing, and equilibration, contact Technical Service.

Void Volume (V_0) Determination:

- Dissolve the Blue Dextran in equilibration buffer containing 5% glycerol at a concentration of 2 mg/mL.
- This concentration of Blue Dextran will give an A_{280} of ~1.0 in the peak fraction.
- Glycerol is added to increase the density of the solution, but its use is optional.
- The recommended sample volume is less than 2% of the total gel bed volume.
- Carefully apply the Blue Dextran sample to the column (avoid disturbing the gel bed surface) to determine V_0 and to check the column packing.
- Immediately after applying the sample, begin collecting fractions of 0.5-1.5% of the total gel bed volume. The flow rate should be ~7% of the column volume per hour.
- Skewing of the Blue Dextran band represents a fault in the column, although some tailing is normal.
- The leading peak indicates the void volume.
- Determine spectrophotometrically the elution volume for Blue Dextran (V_0 for the column) at 280 nm or 610 nm by measuring the volume of effluent collected from the point of sample application to the center of the effluent peak.
- **Note:** Mixing Blue Dextran with kit standards or sample proteins is **not** recommended, since many proteins bind to Blue Dextran.
- Prepare the protein standards and Blue Dextran fresh.²⁸ Occasionally some aggregated protein may appear at the void volume.

Elution Volume (V_e) Determination for Protein Standards:

- Dissolve individual protein standards in equilibration buffer containing 5% glycerol (see **Table 1**).
- If, upon reconstitution, any of the protein solutions contain insoluble material, then filter the protein solution through a 0.45 μ m or 0.2 μ m filter. The loss of protein from this filtration is negligible.
- For a 90 cm × 1.6 cm column, the application of 2.0 mL of individual samples at the recommended concentration (**Table 1**) gives an A_{280} of ~1 in the peak fraction.

Table 1: Recommended Protein Concentrations

Protein	Recommended concentration
Bovine Serum Albumin	10 mg/mL
Apo ferritin	10 mg/mL
Thyroglobulin	8 mg/mL
Alcohol dehydrogenase	5 mg/mL
β -amylase	4 mg/mL
Carbonic anhydrase	3 mg/mL

- Apply the protein standards to the column, using the same sample volume and flow rate as used for the Blue Dextran sample.
- The elution of the standard proteins may be followed by absorbance readings at 280 nm.
- Determine the V_e for the protein standards by measuring the volume of effluent collected from the point of sample application to the center of the effluent peak.

Standard Curve: Plot molecular mass vs. V_e/V_o for each respective protein standard on semi-log paper (see **Figure 1**).

Elution Volume (V_e) Determination for an Unknown Protein:

- Apply the unknown sample to the column at an appropriate concentration using the same sample volume, fraction size, and flow rate as used for the Blue Dextran and the protein standards.
- Determine the V_e of the unknown using the same methods applied to the standards.
- Calculate the V_e/V_o for the unknown and determine its molecular mass from the standard curve.

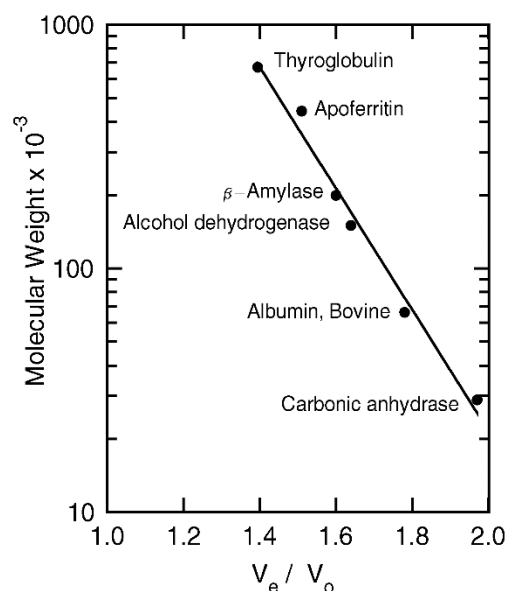


Figure 1. Typical calibration curve obtained with proteins from the MWGF100 Kit run on Sepharose® CL-6B.

Precautions and Disclaimer

For Research Use only. Not for drug, household, or other uses. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

References

1. Whitaker, J.R., *Anal. Chem.*, **35(12)**, 1950-1953 (1963).
2. Andrews, P., *Biochem. J.*, **91(2)**, 222-233 (1964).
3. Flores-Cadengo, Cristina, "Incorporation of EGFR and RON Receptors into Nanodiscs". University of New Mexico, M.S. thesis, p. 20 (May 2019).
4. Ross, Heather A., "Sucrose metabolism in storage organs of *Solanum tuberosum L.*, *Vicia faba L.*, and *Beta vulgaris L.*". University of St. Andrews, Ph.D. dissertation, p. 192, 200 (March 1994).
5. Lipscombe, Richard John, "Mannose Binding Protein Deficiency: Immunochemistry and Mutation Analysis". University of London, Ph.D. dissertation, p. 45 (1995).
6. Arbona, Ricardo Martí, "Mechanistic Characterization of Members of the Amidohydrolase Superfamily". Texas A&M University, Ph.D. dissertation, pp. 24, 88 (December 2006).
7. Lee, Martin Bryce, "Betaine Homocysteine Methyltransferase, Disease and Diet: The Use of Proton Nuclear Magnetic Resonance on Biological Methylamines". University of Canterbury, Ph.D. dissertation, p. 43 (2006).
8. Seyit, Gönül, "Structural Studies of Tripeptidylpeptidase II: Expression and Crystallization Trials". Technische Universität München, Dr. rer. nat. dissertation, p. 41 (2006).
9. Taylor, Alexander I., "Structural and Functional Characterisation of Avian IgY". University of London, Ph.D. dissertation, p. 77 (August 2007).
10. Flodrová, Dana, "Hemicellulose-pectin linkage and exopectate hydrolases involved in homogalacturonan degradation". Brno University of Technology, Ph.D. dissertation, p. 34 (2008).
11. Vanderwerf, Scott, "Signaling Functions of Fanconi Anemia Proteins". Oregon Health Sciences University, Ph.D. dissertation, p. 86 (Summer 2008).
12. Ge, Qiang, "Effect of Lipid Modification and AcrA-Lipid Interactions on the Structure and Oligomerization of AcrA". University of Oklahoma, Ph.D. dissertation, p. 28 (2009).

13. Putz, Annika, "Secondary metabolites from marine sponges, with focus on the chemical ecology and biochemical characterisation of the stress-induced biotransformation of *Aplysina* alkaloids". Heinrich-Heine-Universität Düsseldorf, Ph.D. dissertation, p. 90 (March 2009).
14. Bley, Christopher James, "Mapping the RNA-Protein Interface in Telomerase RNP". Arizona State University, Ph.D. dissertation, p. 61 (May 2011).
15. Chikwana, Vimbai Masiyanise, "Discovery of Novel Amidotransferase Activity Involved In Archaeosine Biosynthesis and Structural and Kinetic Investigation of QueF, an Enzyme Involved in Queuosine Biosynthesis". Portland State University, Ph.D. dissertation, p. 23 (2011).
16. Nicholls, Samantha Elizabeth Bernard, "Development and Characterization of Caspase Activatable GFP and a Family of Fluorescent Reporters". University of Massachusetts Amherst, Ph.D. dissertation, p. 71 (February 2013).
17. Wetter, Michael Lukas, "Development of Novel Conjugate Vaccines Against *Salmonella*". Universität Zürich, Dr. sc. nat. dissertation, p. 145 (April 2014).
18. Sun, Tao, "*Trans*-acting factors essential for plant organelle RNA editing". Cornell University, Ph.D. dissertation, p. 71 (2014).
19. Greene, Landon Ray, "Isolation, Identification, and Quantitation of Biomolecules using High Performance Liquid Chromatography and Mass Spectrometry". Rutgers, The State University of New Jersey, Ph.D. dissertation, p. 92 (October 2014).
20. Zhao, Yuting, "The Role of the Exocyst in Membrane Deformation, Cell Migration and Exocytosis and Exocytosis". University of Pennsylvania, Ph.D. dissertation, p. 66 (2014).
21. Tekin, Izel, "Genetic contributors to non-motor symptoms of sporadic Parkinson's Disease". The Pennsylvania State University, Ph.D. dissertation, p. 78 (May 2016).
22. Petersburg, Jacob Richard, "Prosthetic antigen receptors as a platform for solid tumor immunotherapy". University of Minnesota, Ph.D. dissertation, pp. 199, 221 (December 2017).
23. Liu, Feng, "Functional characterization of mitochondrial and cytosolic aldehyde dehydrogenases in maize (*Zea mays L*)". Iowa State University, Ph.D. dissertation, pp. 136, 198 (2017).

24. Liu, Yi, "Characterization of Extracellular Polymeric Substances in Biological Wastewater Treatment Systems". University of Guelph, Ph.D. dissertation, p. 36 (May 2018).
25. Weijman, Johannes F., "Biochemistry of Apoptosis Signal-regulating Kinase 1 Regulation and Hydrogen Peroxide Signalling". University of Otago, p. 213 (October 2018).
26. Rai, Shashank, "Role of WD domain of ATG16L1 in maintaining autophagy and tissue homeostasis in mice". University of East Anglia, Ph.D. dissertation, p. 43 (2019).
27. McWhite, Claire Darnell, "Conservation and comparison of protein interactions across evolution". University of Texas at Austin, Ph.D. dissertation, p. 82 (May 2020).
28. Marshall, J.J., *J. Chromatog.*, **53(2)**, 379-380 (1970).

Notice

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

Technical Assistance

Visit the tech service page at [SigmaAldrich.com/techservice](https://www.sigmaaldrich.com/techservice).

Terms and Conditions of Sale

Warranty, use restrictions, and other conditions of sale may be found at [SigmaAldrich.com/terms](https://www.sigmaaldrich.com/terms).

Contact Information

For the location of the office nearest you, go to [SigmaAldrich.com/offices](https://www.sigmaaldrich.com/offices).

The life science business of Merck KGaA, Darmstadt, Germany operates as MilliporeSigma in the U.S. and Canada.

MilliporeSigma and Sigma-Aldrich are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.

© 2015-2026 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved.

23357251 Rev 03/26

**MILLIPORE
SIGMA**