

Version 1a Last updated 27 July 2021

ab193259 Protein G Sepharose®

For the purification of monoclonal and polyclonal antibodies.

View Protein G Sepharose® Column datasheet:

www.abcam.com/ab193259

[use www.abcam.cn/ab193259 for China, or www.abcam.co.jp/ab193259 for Japan]

This product is for research use only and is not intended for diagnostic use.

Table of Contents

1. Overview	1
2. Materials Supplied and Storage	2
3. General guidelines, precautions, and troubleshooting	2
4. Assay Procedure	3
5. Notes	4

1. Overview

Protein G is a cell wall protein produced by group G streptococcus. Like protein A, this bacteria-derived protein binds with high affinity and specificity to the Fc portion of most mammalian immunoglobulins. Therefore, Protein G has been widely used for IgG purification.

Protein G Sepharose® ab193259 is a genetically engineered protein containing three Ig-binding regions of native Protein G. The cell wall binding region, albumin binding region and other non-specific regions have been eliminated from the recombinant Protein G to ensure the maximum specific IgG binding. The coupling technique is optimized to give a higher binding capacity for IgG and minimum leaching of recombinant Protein G. In addition, Protein G-Sepharose beads display high chemical and physical stability as well as high flow rate, hydrophilicity and high gel strength. It can be used for IgG purification and immunoprecipitation. 6% cross-linked Sepharose beads supplied as 50% slurry (e.g., 1 ml of settled resin is equivalent to 2 ml of 50% slurry) in 20% Ethanol/H₂O

Flow rate tested: 2.07 mL/min

Test condition: Calculations based on the time required to pass 18 ml of water through 2 ml settled beads (column diameter 1.5 cm).

Usage: Reusable for up to 10 times without significant loss of binding capacity.

2. Materials Supplied and Storage

Store at 4°C. Do not freeze. Stable, as supplied, for at least 1 year.

3. Materials Required but not supplied

Binding Buffers: PBS/TBS/0.15 M sodium chloride in 50 mM sodium borate, pH 8.0.

Elution Buffers: 0.1 M citric acid, pH 2.75.

Neutralization Buffer: 1 M Tris-HCl, pH 9.

Ready-to-use pre-packed columns.

4. General guidelines, precautions, and troubleshooting

Please observe safe laboratory practice and consult the safety datasheet.

For general guidelines, precautions, limitations on the use of our assay kits and general assay troubleshooting tips, particularly for first time users, please consult our guide:

www.abcam.com/assaykitguidelines

For typical data produced using the assay, please see the assay kit datasheet on our website.

5. Assay Procedure

- 1.1 Carefully pack the column avoiding air bubbles.
- 1.2 Equilibrate the column with 5X resin bed volume of Binding Buffer and allow the buffer to drain through the column. Do not let the resin bed dry.
- 1.3 Dilute serum sample with Binding Buffer (1:1 ratio).
- 1.4 Mix the diluted serum sample well. Make sure there are no bubbles in the sample solution.
- 1.5 Apply the diluted sample onto the column. Do not let the resin bed dry.
- 1.6 Collect the flow-through.
- 1.7 Reapply the flow-through to the column and collect the sample. Repeat 4 times.
- 1.8 Wash the column 4 – 5 times with 5X volume of Binding Buffer containing 0.5 M NaCl.
- 1.9 Wash the column 4 - 5 times with Binding Buffer.
- 1.10 Elute antibodies with Elution Buffer ~3-5X resin bed volume.
- 1.11 Collect fractions using micro centrifuge tube. Immediately neutralize the eluted fractions by adding 100 μ L of 1 M Tris, pH 9.0 per mL of eluate.
- 1.12 Assay protein concentration by measuring the absorbance at 280 nm and combine the fractions with highest absorbance.
1 OD₂₈₀ = 0.73 mg/mL IgG.
- 1.13 To regenerate/store column:
 - 1.13.1 Wash with 5 volumes of Elution Buffer.
 - 1.13.2 Wash with 5 volumes of distilled water.
 - 1.13.3 Store column in 20 % Ethanol/H₂O at 4°C. Store upright at 4°C.

6. Notes

Technical Support

The Abcam logo is a registered trademark.
All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)