

## ab193262 – Protein A/G Sepharose®

For the purification of monoclonal and polyclonal antibodies.

For overview, typical data and additional information please visit:

<http://www.abcam.com/ab193262>

### Procedure Example

- 1) Wash column with ddH<sub>2</sub>O to remove air bubbles.
- 2) Fill column with protein A/G beads.
- 3) Wash the column with 5X volume of Binding Buffer.
- 4) Dilute serum sample with Binding Buffer (1:1 ratio).
- 5) Invert the diluted serum sample to mix well. Make sure no bubbles in the solution.
- 6) Pour the solution onto the column.
- 7) Collect the solution and repeat step 6 & 7 for 10 times.
- 8) Wash the column 4 – 5 times with Binding Buffer containing 0.5 M NaCl.
- 9) Wash the column 4 - 5 times with the Binding Buffer.
- 10) Add Elution Buffer to elute IgG (0.5-1 ml each time).
- 11) Collect the eluent using microcentrifuge tube.
- 12) Assay protein concentration and combine the fractions containing sufficient amount of IgG.
- 13) To regenerate/store column:
  - a. Wash with 3 volumes of elution buffer.
  - b. Wash with 3 volumes of distilled water.
  - c. Store column in 20 % Ethanol/H<sub>2</sub>O.

### Buffer Example

- 1) Binding buffer: 0.05 M sodium borate, 0.15 M sodium chloride pH 8.0.
- 2) Elution buffer: 0.1 M citric acid, pH 2.75.

### Technical Support

Copyright © 2023 Abcam. All Rights Reserved. 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](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)