

## ab288103 – Protein A Antibody Purification Kit

A simple, ready to use kit containing all necessary buffers and pre-packed columns to purify up to 350 mg of IgG.  
For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit:

<http://www.abcam.com/ab288103>

### Storage and Stability

Store kit at 4°C, protected from light.

### Materials Supplied

Item	Quantity	Storage Condition
Hi-Bind™ Protein Column	1 column	+4°C
Binding Buffer (10X)	40 ml	+4°C
Elution Buffer (10X)	20 ml	+4°C
Neutralization Buffer	10 ml	+4°C

### Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully utilize this assay:

- 20% Ethanol
- 1.5 ml Collection Tubes
- 50 ml Centrifuge Tubes

### Reagent Preparation

- Briefly centrifuge small vials prior to opening.
- Read the entire protocol before purifying IgG.

**Binding Buffer:** Dilute Binding Buffer (10X) with ddH<sub>2</sub>O. (i.e. add 1 ml Binding buffer (10X) to 9 ml ddH<sub>2</sub>O). Diluted buffer is stable at 4°C for 3 months.

**Elution Buffer:** Dilute Elution Buffer (10X) with ddH<sub>2</sub>O. (i.e. add 1 ml Elution buffer (10X) to 9 ml ddH<sub>2</sub>O). Diluted buffer is stable at 4°C for 3 months.

### Antibody Purification Protocol

1. Centrifuge samples at 10000 x g at 4°C for 25 min and collect the supernatant to a new tube. Equilibrate samples by mixing 1:1 (v/v) of sample with Binding Buffer. The maximum loading volume is 4 ml at once. (For larger sample size, equilibrate samples by mixing 9 volume of sample with 1 volume of 10X Binding Buffer.)
2. Remove the top cap and bottom cap respectively. Allow the column to drain the storage buffer. Equilibrate the column by adding 5 ml Binding Buffer. Do not let resin bead dry out at any time.
3. Place a clean 50 ml Centrifuge Tube underneath the Column. Gradually load the equilibrated sample into the Column and let it flow through the Column. Reload the non-adsorbed material 8 to 10 times at RT to maximize the IgG binding capacity.
4. Collect the final flow through. Wash the Column twice with 5 ml Binding Buffer.

**ΔNote:** *Keep the flow through until satisfactory enrichment of IgG in eluate is confirmed.*

5. Prepare 10 collection tubes (label 1-10) with 100 µl Neutralization Buffer in each tube. Put the Column in tube #1 and add 1 ml Elution Buffer in the Column. Mix the eluted

samples with the Neutralization Buffer immediately. Repeat the elution step for another 6-9 times, each time with a new Collection Tube.

6. **Measurement:** Measure the IgG concentration for each fraction by measuring absorbance at 280 nm (1.4 OD<sub>280</sub> ≈ 1 mg/ml IgG). Combine the eluted fractions containing purified IgG.
7. **Column Regeneration:** Wash the Column with 10 ml Elution Buffer, followed by 5 ml distilled water twice and then 5 ml Binding Buffer twice. Add 5 ml 20% ethanol into the Column, drain out ~3 ml, and store the Column at 4 °C. Column can be regenerated up to 10 times without significant loss of binding capacity. We recommend reusing the column only for the same sample to avoid cross contamination.

### Technical Support

Copyright © 2021 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)