

ab286842 – Protein A/G Magnetic Beads

For Purification of antibodies from multiple sources.
For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit:

<http://www.abcam.com/ab286842>

Storage and Stability

On receipt product should be stored at +4°C. Do not freeze. Upon opening, use kit within 12 months.

Materials Supplied

Protein A/G Magnetic Beads supplied as a 50% slurry in PBS with 0.02% sodium azide.

Technical Specifications

Support Characteristics: Paramagnetic, spherical, 6% cross-linked agarose.

Ligand: Recombinant fusion Protein A/G.

Particle size: 75 – 150 µm.

Binding capacity: Generally >10 mg human IgG/ml wet beads.

Procedure Protocol

Δ Note: Prepare the antibody solution by diluting the required amount of antibody in binding buffer before running the protocol.

Magnetic bead preparation:

1. Dispense the required amount of magnetic beads into 1.5 ml microfuge tube.
2. Place tube in magnetic rack and remove storage solution.
3. Add 500µl binding buffer.
4. Resuspend the beads.
5. Remove liquid.

Antibody capture:

1. Immediately add the antibody solution.
2. Resuspend and mix (slow end-over-end) for at least 15 minutes.
3. Remove the liquid.

Washing:

1. Add 500 µl Binding Buffer containing 0.5 M NaCl; Remove the liquid.
2. Add 500 µl Binding Buffer; Remove liquid.

Target Binding:

1. Add sample diluted in binding buffer.
2. Incubate with slow end-over-end mixing for up to 60 minutes.
3. Remove and collect unbound fraction.

Washing: (perform 3 times)

1. Add 500 µl wash buffer.
2. Remove liquid (save washes to troubleshoot).

Elution: (perform 3 times)

1. Add 2 volumes elution buffer (vs. bead volume).
2. Completely resuspend beads and incubate for at least 2 minutes.
3. Remove and collect elution fraction.

Buffer examples:

(1) Binding Buffer: 50 mM Tris, 150mM NaCl, pH 7.5.

(2) Wash Buffer: 50 mM Tris, 150mM NaCl, pH 7.5 (Or add 1% Octyl glucoside to this buffer)

(3) Elution Buffer: 0.1 – 0.2M Glycine pH 2.5-3.1 (or 0.1M citric acid, pH 2.5-3.1 or 2.5% Acetic Acid)

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

www.abcam.cn/contactus (China)

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