

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:

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 20% slurry in PBS with 0.02% sodium azide.

Technical Specifications

Ligand: Recombinant fusion Protein A/G.

Particle size: 30 – 100 µm.

Binding capacity: 10 mg Rabbit IgG/ml beads

Procedure Protocol

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

Δ Note: This protocol is designed for 100 µL Protein A/G Magnetic Beads but it may be adjusted as required.

Magnetic bead preparation:

1. Resuspend the beads by shaking or vortexing the vial.
2. Transfer 500 µl Protein A/G Magnetic Beads into a clean tube.
3. Place the tube on a magnetic separation rack to collect the beads. Remove and discard the supernatant.
4. Add 1 ml Binding/Wash Buffer to the tube and invert the tube several times to mix. Use the magnetic separation rack to collect the beads and discard the supernatant. Repeat this step twice.

Protein purification:

1. Resuspend the beads in 100 µl Binding/Wash Buffer.
2. Add the sample to the tube and gently invert the tube to mix.
3. Incubate the tube at room temperature with mixing (on a shaker or rotator) for 30 minutes. It also can be done at 4°C overnight.
4. Use the magnetic separation rack to collect the beads and discard the supernatant. If necessary, keep the supernatant for analysis.
5. Add 1ml Binding/Wash Buffer to the tube and mix well, use the magnetic separation rack to collect the beads and discard the supernatant. Repeat the wash step three more times.

Elution: (perform 3 times)

1. Add 300-500 µL elution buffer, mix well and incubate for 5-10 minutes on a shaker or rotator.
2. Use the magnetic separation rack to collect the beads and transfer the supernatant into a new tube. Repeat this step.
3. Add 50 µl Neutralization buffer per 500 µl eluate to neutralize the pH.

Buffer examples:

Binding/Wash Buffer: PBS, pH 7.4

Elution Buffer: 0.1M Glycine pH 2-3

Neutralization buffer: 1M Tris, pH 8.5

For technical support contact information, visit: www.abcam.com/contactus