

ab286896 – Endotoxin-Free Protein A Agarose

For purification of monoclonal and polyclonal antibodies from culture media, serum, ascites fluid or hybridoma supernatants.
For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit:

<http://www.abcam.com/ab286896>

Protocol Example (Antibody Purification)

1. Carefully pack the column avoiding air bubbles.
2. Equilibrate the column with 5 resin volume of endotoxin-free Binding Buffer (i.e. PBS) & allow the buffer to drain through the column. Do not let the resin dry.
3. Dilute serum sample with endotoxin-free Binding Buffer (1:1 ratio).
4. Mix well the diluted serum sample. Make sure there are no bubbles in the sample solution.
5. Apply the diluted sample onto the column. Do not let the resin dry.
6. Collect the flow-through.
7. Reapply the flow-through to the column & collect the sample. Repeat 4 times.
8. Wash the column 4 - 5 times with 5 volume of endotoxin-free Binding Buffer containing 0.5 M NaCl.
9. Wash the column 4 - 5 times with endotoxin-free Binding Buffer.
10. Elute antibodies with endotoxin-free Elution Buffer (i.e. 0.1 M Citric Acid) ~3-5 resin volume. Collect fractions using micro centrifuge tube containing endotoxin-free neutralization buffer (100 μ l of 1 M Tris, pH 9.0 per ml of eluate).
11. Assay protein concentration by measuring the absorbance at 280 nm and combine the fractions with highest absorbance. 1 OD₂₈₀ = 0.73 mg/ml IgG.

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)