

ab128745 Mouse Antibody Purification Kit

A product of Expedeon, an
Abcam company

Applicable to Expedeon product codes 830-0005, 830-0010.

View ab128745

Mouse Antibody Purification Kit datasheet:

www.abcam.com/ab128745

(use www.abcam.cn/ab128745 for China, or www.abcam.co.jp/ab128745 for Japan)

For preparing antibodies for conjugation.

This product is for research use only and is not intended for
diagnostic use.

Table of Contents

| | |
|-----------------------------------|---|
| 1. Overview | 2 |
| 2. Materials Supplied and Storage | 3 |
| 3. Technical Considerations | 4 |
| 4. Assay Procedure | 5 |

1. Overview

Commercially available antibodies often contain substances (e.g. BSA, glycine, Tris, azide) that interfere in labeling reactions with enzymes or fluorophores. ab128745 quickly removes these contaminants. It can also be used to purify antibodies from crude samples such as ascites fluid or immune serum.

The resin has a high affinity for mouse IgG molecules. The method involves capture of the antibody on a Mouse resin and the removal of unwanted substances by a simple wash procedure, which is carried out in a standard microfuge. The purified product is then eluted and neutralized.

Δ Note: *The Mouse Antibody Purification Kit is not suitable for use with antibodies from other species. It should be noted that the binding strength of bovine IgG to the Mouse resin is negligible, and therefore this kit is especially suitable for purification of antibodies from cell culture that has contained bovine calf serum.*

Antibodies purified using the Mouse Antibody Purification Kit are fully compatible with our [Lightning-Link® Antibody Conjugation kits and our Oligonucleotide Conjugation Kit](#).

2. Materials Supplied and Storage

Store at +4°C upon receipt. Do not freeze or store the resin at room temperature.

| Item | Quantity | | Storage temperature |
|---|----------|----------|---------------------|
| | 1 x Test | 3 x Test | |
| Spin Cartridge/Collecting Tube Assembly | 1 unit | 3 units | +4°C |
| Additional Collecting Tubes | 4 units | 12 units | +4°C |
| 10x Binding Buffer | 1 vial | 1 vial | +4°C |
| Wash Buffer | 1 vial | 1 vial | +4°C |
| Elution Buffer | 1 vial | 1 vial | +4°C |
| Mouse resin | 1 vial | 3 vials | +4°C |
| Neutralization Buffer | 1 vial | 1 vial | +4°C |

Reagents are ready to use as supplied.

3. Technical Considerations

Recommended antibody quantities:

The antibody to be purified or cleaned up is ideally in a volume of 100 μ L to 0.5 mL. Up to 150 μ g of antibody can be purified in each run.

Antibody pre-conjugation considerations:

This kit can be used for preparing antibodies for conjugation. The antibody concentration for each Conjugation Kit has been optimised. Before starting the elution step of this purification procedure, please refer to the relevant Lightning-Link® Conjugation Kit datasheet or protocol for the recommended antibody concentration and find more general information about antibody conjugation at www.abcam.com/conjugationFAQs.

Test for protein concentration:

Wherever possible, protein values should be determined using an absorbance at 280 nm. An extinction co-efficient of 1.4 is generally used for IgG – so a 1 mg/mL solution of IgG will give an absorbance value of 1.4 when measured with a 1 cm path length.

Δ Note: *If a low volume/amount of antibody has been added, the concentration of protein in the eluates will be low.*

When other methods of determining IgG concentration are used such as BCA or Bradford protein assays, determinations should be performed before the addition of the Neutralization Buffer, as this can interfere with these reagents. Remove an aliquot for protein determination and neutralize the rest of the fraction immediately as the low pH of the Elution Buffer can denature the antibody.

When using Bradford-type reagents it is important to use an IgG standard curve. Failure to do this will result in incorrect antibody levels being calculated. If IgG is not available then a BSA standard curve can be used, but the IgG levels will be under-estimated by a factor of 2.3.

4. Assay Procedure

Reconstitution of Mouse Resin:

Add 0.3 mL of Wash Buffer to each vial of Mouse Resin, mix by inversion for a few seconds and transfer to the spin cartridge. Spin for 30 seconds in a microfuge.

Incubation of Sample with Resin:

To the antibody, add an appropriate amount of 10x Binding Buffer which corresponds to 1/10th of the sample volume. For example, if the sample volume is 200 μ L, add 20 μ L of Binding Buffer. Pipette the sample into the spin cartridge and cap the tube. Incubate for a minimum of 2 hours with agitation or periodic shaking. Alternatively, incubate overnight at either +4°C or room temperature.

***Δ Note:** The volume of antibody to be purified or cleaned up should ideally be 0.1-0.5 mL, though larger volumes may be processed by first incubating the antibody sample (combined with Binding Buffer) with the Mouse resin in a larger vessel (e.g. 2 mL tube) prior to transferring to the spin cartridge in several aliquots, spinning own excess liquid each time.*

For larger volumes (cell culture supernatant) reconstitute the resin with 0.3 mL of wash buffer, transfer to the spin cartridge and centrifuge to get rid of the liquid (this can be discarded). Add 1/10 volume of binding buffer (BB) to the cell supernatant containing the antibody of interest. Add ~0.4 mL of supernatant + BB to the resin, mixing well and transferring the whole volume into the remaining sample. This process can be done twice to make sure all the resin is collected from the spin cartridge.

Wash Procedure:

Microfuge the spin cartridge assembly for 30 seconds to remove most of the non-bound protein. Add 0.5 mL of Wash Buffer and spin again. Repeat the wash procedure three times.

***Δ Note:** Save the non-bound and wash fractions by transferring the material from the collecting tube after each spin to a set of tubes (not supplied). Do not use the four*

collecting tubes supplied with the kit, as these have an extended hinge to accommodate the spin cartridge, and are required for the elution step.

Elution:

Please see the Technical Considerations sections 3.2 and 3.3 before starting this step.

Transfer the cartridge to a clean collecting tube. Add 100 μ L of Elution Buffer and incubate for 2 minutes at room temperature with gentle agitation. Microfuge for 30 seconds. Remove the collecting tube and add 25 μ L Neutralization Buffer to the tube.

Place the cartridge in a new collecting tube and add a further 100 μ L of Elution Buffer to the Mouse resin. Incubate for 2 minutes at room temperature with gentle agitation. Spin and collect and neutralize as before.

Repeat the elution procedure until all four clean collecting tubes have been used. The protein normally elutes in tubes 1 and 2 but you should confirm this using a test for protein concentration (see Technical Consideration section 3.3) before pooling any of the tubes.

Pool the tubes with most protein (normally two tubes; if more than two tubes are strongly positive it is possible that you have used too much sample in your protein assay). However, if your application does not require a high concentration of antibody you may choose to pool all tubes that contain protein, regardless of concentration.

Antibody storage:

Store at +4°C. Other storage conditions (e.g. frozen at -70°C) may also be satisfactory). The sensitivity of any particular mouse antibody to freeze-thaw should be determined by experimentation on small aliquots.

Version 6 Last Updated February 5, 2024

Technical Support

Copyright © 2024 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)