

Antibody Purification Kit (Protein A) ab102784

4 References

Overview

Product name Antibody Purification Kit (Protein A)

Product overview Commercially available antibodies often contain substances (e.g. BSA, glycine, tris, azide) that interfere in labeling reactions with enzymes or fluorophores. **ab102784** quickly removes these contaminants. It can also be used to purify antibodies from crude samples such as ascites fluid or immune serum. The antibody to be purified or cleaned up is ideally in a volume of 100-500 µl.

Up to 20-500 µg of antibody can be purified in each run.

The method involves capture of the antibody on Protein A 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.

Note: This method cannot be used with samples containing relatively dilute antibody in large volumes (e.g. tissue culture supernatant). For larger volumes, we recommend the use of Antibody TCS Purification Kit (3 purifications) ([ab109207](#)) or Antibody Serum Purification kit (3 purifications) ([ab109209](#)).

Antibodies purified using the Antibody Purification Kit (Protein A) are fully compatible with our Lightning-Link® Antibody Conjugation Kits and our Oligonucleotide Conjugation Kit. However they are not compatible with our GOLD Antibody conjugation kits. To purify antibodies for use with our Gold conjugation kits, please use our Gold antibody purification kit ([ab204909](#))

ab102784 is not suitable for goat antibody purification. Please refer to the protocol booklet for more information about suitable and unsuitable species antibodies."

Notes This product is manufactured by Expeleon, an Abcam company, and was previously called AbSelect™ Antibody Purification System. 860-0005 is the same as the 1 test size. 860-0010 is the same as the 3 tests size.

Properties

Storage instructions Store at +4°C. Please refer to protocols.

Components	3 tests	1 tests
ab273959 - Additional Collecting Tubes	12 vials	4 vials
Binding Buffer (10X)	1 x 200µl	1 x 200µl
Elution Buffer	1 x 1500µl	1 x 1500µl
Neutralizer Buffer	1 x 500µl	1 x 500µl
Protein A resin	3 x 100µl	1 x 100µl
ab273958 - Spin cartridge/ collecting tube assemblies	3 units	1 unit
Wash Buffer	1 x 12ml	1 x 12ml

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors