

Antibody Purification Kit (Protein G) ab128747

1 References

概述

产品名称	抗体Purification试剂盒(Protein G)
产品概述	<p>Commercially available antibodies often contain substances (e.g. BSA, glycine, tris, azide) that interfere in labeling reactions with enzymes or fluorophores. ab128747 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 should be in a volume of 100 to 500 µL. 20 to 300 µg of antibody can be purified in each run.</p> <p>The method involves capturing the antibody on the Protein G resin. Protein G has a high affinity for the Fc regions of IgG molecules from a variety of species. Once the antibody has bound to the Protein G, unwanted substances can be removed by simply washing the resin. Up to 300µg of antibody can be purified in each run. The purified product is then eluted and neutralized.</p> <p>Antibodies purified with the Antibody Purification Kit (Protein G) are fully compatible with our Lightning-Link® Antibody Conjugation kits and Oligonucleotide Conjugation Kit. To purify antibodies for use with our Gold conjugation kits, please use our Gold antibody purification kit (<a href="#">ab204909</a>).</p>
说明	<p>This product is manufactured by Expedeon, an Abcam company, and was previously called AbSelect™ G Antibody Purification System. 890-0005 is the same as the 1 test size. 890-0010 is the same as the 3 tests size.</p>

性能

存放说明	Store at +4°C. Please refer to protocols.
------	---

组件	1 tests	3 tests
ab273959 - Additional Collecting Tubes	4 vials	12 vials
Elution Buffer	1 x 1500µl	1 x 1500µl
ab275847 - G Binding Buffer (10X)	1 x 200µl	1 x 200µl

组件	1 tests	3 tests
Neutralizer Buffer	1 x 500µl	1 x 500µl
ab275848 - Protein G Resin	1 x 400µl	3 x 400µl
ab273958 - Spin cartridge/ collecting tube assemblies	1 unit	3 units
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.cn/abpromise> or contact our technical team.

### Terms and conditions

---

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