

Serum Antibody Purification Kit (Protein G) ab128751

概述

产品名称 Serum抗体Purification试剂盒(Protein G)

样品类型 Serum, Ascites Fluid

产品概述 Protein G has a high affinity for the Fc regions of IgG molecules from a variety of species. ab128751 G Serum resin is prepared by coupling purified Protein G to agarose beads. It can therefore be used to purify IgG fractions from both serum and ascites fluid. The antibody is captured on the AbSelect G resin and unwanted substances are removed by a simple wash procedure. Up to 10mg of antibody can be purified in each run. The volume of sample required will depend on the host species.

The purified product is then eluted and neutralized.

The antibodies purified with Protein G Serum Antibody Purification Kit are fully compatible with our Lightning-Link[®] antibody conjugation kits and Oligonucleotide Conjugation Kit. To purify antibodies for use with our nanoparticle conjugation kits, please use Antibody Purification Kit - Nanoparticles ([ab204909](#)).

说明 This product is manufactured by Expedeon, an Abcam company, and was previously called AbSelect™ G Serum Antibody Purification System. 893-0030 is the same as the 1 test size. 893-0500 is the same as the 3 tests size.

性能

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

组件	1 tests	3 tests
Elution Buffer	1 x 5ml	1 x 15ml
ab275847 - G Binding Buffer (10X)	1 x 6ml	1 x 18ml
Neutralizer Buffer	1 x 2ml	1 x 5ml
ab274180 - Purification Column	1 unit	3 units

组件	1 tests	3 tests
ab275849 - Serum Protein G Resin	1 x 4ml	3 x 4ml
Wash Buffer	1 x 25ml	1 x 75ml

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