

Serum Antibody Purification Kit (Protein G) ab128751

製品の概要

製品名	Serum Antibody Purification Kit (Protein G)
サンプルの種類	Serum, Ascites Fluid
製品の概要	<p>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.</p> <p>The purified product is then eluted and neutralized.</p> <p>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).</p>
特記事項	<p>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.</p>

製品の特性

保存方法 Store at +4°C. Please refer to protocols.

内容	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

内容	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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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