

TCS Antibody Purification Kit ab109207

概述

产品名称

TCS抗体Purification试剂盒

产品概述

TCS Antibody Purification Kit (ab109207) has been designed for tissue culture supernatant purification, and can be used to purify 10-50 mL in each purification. Each column can purify up to 5 mg of antibody.

This kit works by coupling highly purified protein A to agarose beads and can therefore be used to purify IgG fractions from hybridoma supernatants. The antibody is captured on the TCS resin and unwanted substances are removed by a simple wash procedure. The purified product is then eluted and neutralized.

The antibodies purified with the TCS Antibody Purification Kit are fully compatible with our Lightnign-Link[®] Antibody Conjugation kits and our Oligonucleotide Conjugation Kit. To purify antibodies from TCS for use with our nanoparticle conjugation kits, please use our TCS Antibody Purification Kit - Nanoparticles ([ab270554](#)).

ab109207 is not suitable for goat antibody purification.

说明

This product is manufactured by Expedeon, an Abcam company, and was previously called AbSelect[™] A Antibody Purification System. 862-0500 is the same as the 3 tests size. 862-0030 is the same as the 1 test size.

性能

存放说明

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

组件	3 tests	1 tests
Binding Buffer (10X)	1 x 18ml	1 x 6ml
Elution Buffer	1 x 15ml	1 x 5ml
Neutralizer Buffer	1 x 5ml	1 x 2ml
Protein A resin	3 x 2ml	1 x 2ml

组件	3 tests	1 tests
ab274180 - Purification Column	3 units	1 unit
ab273958 - Spin cartridge/ collecting tube assemblies	3 units	1 unit
Wash Buffer	1 x 75ml	1 x 25ml

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