

TCS Antibody Purification Kit ab109207

Overview

Product name TCS Antibody Purification Kit

Product overview 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 Lightning-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.

Notes 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.

Properties

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

Components	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

Components	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.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