

Mouse TCS Antibody Purification Kit ab128749

[5 References](#) [2 Images](#)

Overview

Product name	Mouse TCS Antibody Purification Kit
Product overview	<p>Efficient mouse antibody purification from tissue culture supernatants. Processes 10-25 mL of supernatant to purify up to 1.5 mg of antibody per run.</p> <p>The mouse resin has a very high affinity and specificity for mouse IgG molecules. ab128749 Mouse TCS Antibody Purification Kit can be used to purify mouse IgG fractions from hybridoma supernatants. The binding strength of bovine IgG to Abcam's Mouse resin is negligible; therefore, it can be used to selectively recover mouse IgG from TCS samples.</p> <p>The method involves capture of the mouse antibody on the Mouse resin and removal of unwanted substances using a simple wash procedure. The purified product is eluted and neutralized.</p> <p>Antibodies purified using the Mouse TCS Antibody Purification kit are fully compatible with our Lightning-Link® Antibody Conjugation Kits and our Oligonucleotide Conjugation Kit, however they are not compatible with our nanoparticle Antibody conjugation kits. To purify antibodies for use with our nanoparticle conjugation kits (Gold, Latex, Europium and Magnetic), please use our <u>Mouse Antibody Purification Kit - Nanoparticles ab204910</u>.</p> <p>ab128749 is suitable for use with Mouse antibodies. It is not suitable for use with antibodies from other species.</p>
Notes	<p>This product is manufactured by Expedeon, an Abcam company, and was previously called AbSelect™ Mouse TCS Antibody Purification System. 832-0500 is the same as the 3 tests size. 832-0005 is the same as the 1 test size.</p>

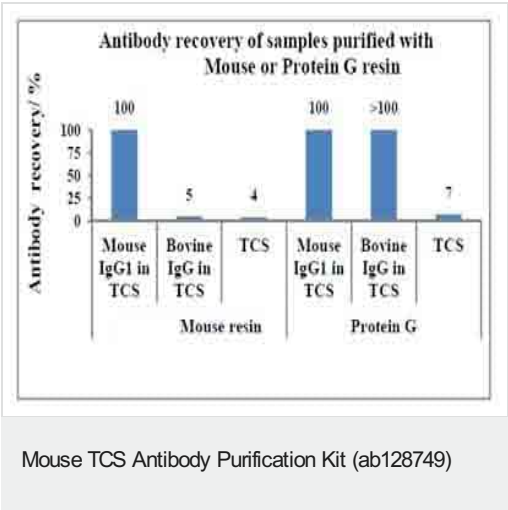
Properties

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

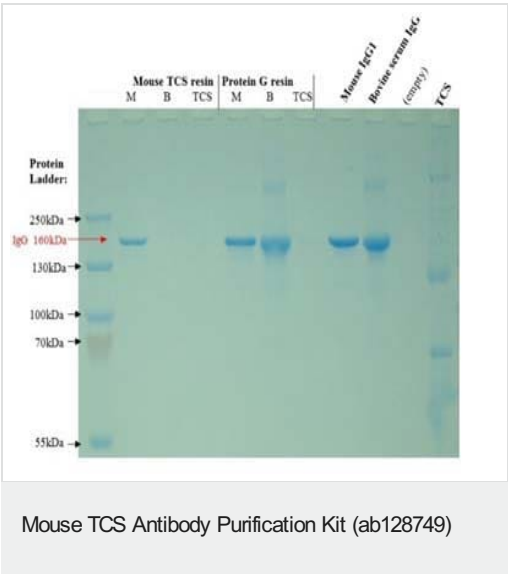
Components	1 tests	3 tests
Binding Buffer (10X)	1 x 6ml	1 x 18ml

Components	1 tests	3 tests
Elution Buffer	1 x 5ml	1 x 15ml
ab273960 - Mouse Resin	1 x 2ml	3 x 2ml
Neutralizer Buffer	1 x 2ml	1 x 5ml
ab274180 - Purification Column	1 unit	3 units
ab273914 - Spin Cartridge / Collecting Tube Assemblies	1 unit	3 units
Wash Buffer	1 x 25ml	1 x 75ml

Images



A representative bar graph comparing the effectiveness of Mouse TCS Antibody Purification Kit and Protein G resin.



Gel electrophoresis of TCS samples purified with the Mouse TCS Antibody Purification Kit.

- M = Mouse IgG1 in a TCS sample
- B = Bovine serum IgG in a TCS sample
- TCS = Tissue culture supernatant sample

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