

Product datasheet

Rat TCS Antibody Purification Kit ab270562

1 Image

Overview

Product name Rat TCS Antibody Purification Kit

Product overview Rat TCS Antibody Purification Kit (ab270562) is designed for the purification of rat IgG fractions from tissue culture supernatants. The Rat Resin has a very high affinity and specificity for rat IgG molecules. The binding strength of bovine IgG to the Rat Resin is negligible; therefore, it can be used to selectively recover rat IgG from TCS samples. The rat antibody to be purified should be in 10 to 50 mL of tissue culture supernatant. Up to 600 µg of antibody can be purified in each run.

This product is manufactured by Expedeon, an Abcam company. It was previously called AbSelect™ Rat TCS Antibody Purification System. Expedeon product code 842-0500 is the same as the 3 pack size of this product.

The method involves capture of the rat antibody on the Rat resin and removal of unwanted substances using a simple wash procedure. The purified product is eluted and neutralized. Rat TCS Antibody Purification Kit is not suitable for use with antibodies from other species.

Antibodies purified using the Rat TCS Antibody Purification Kit are fully compatible with Lightning-Link® Antibody Conjugation kits and our Oligonucleotide Conjugation Kit.

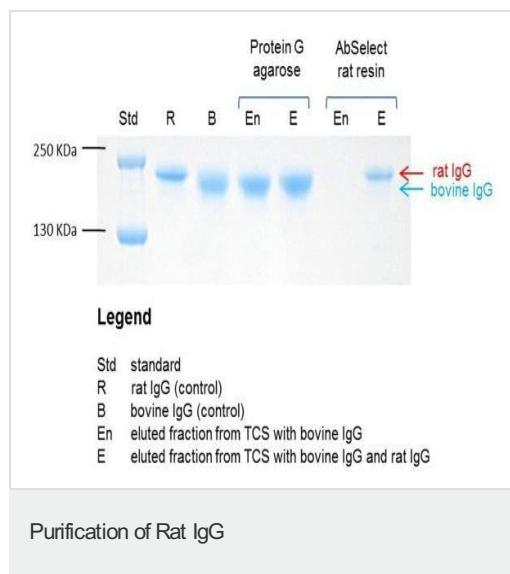
Properties

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

Components	3 tests
ab274185 - AbSelect	3 x 400µl
ab273959 - Additional Collecting Tubes	12 vials
ab274298 - Binding Buffer (10X)	1 x 18ml
ab274299 - Elution Buffer	1 x 1500µl
ab274174 - Neutralizer Buffer	1 x 500µl

Components	3 tests
ab273914 - Spin Cartridge / Collecting Tube Assemblies	3 units
ab273958 - Spin cartridge/ collecting tube assemblies	3 units
ab274173 - Wash Buffer	1 x 12ml

Images



Purification of rat IgG using ab270562.

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