

TCS Antibody Purification Kit (Gold/Europium labelling) ab270554

画像数 1

製品の概要

製品名 TCS Antibody Purification Kit (Gold/Europium labelling)

製品の概要 TCS Antibody Purification Kit - Nanoparticles (ab270554) is designed for the purification of IgG fractions from tissue culture supernatants. TCS Protein A resin is prepared by coupling purified Protein A to agarose beads.

The method involves capture of the antibody on the TCS Protein A resin and the removal of unwanted substances using a simple wash procedure. The antibody is then eluted and neutralized. The kit can be used in its rapid version by spin-purification in a centrifuge, or by gravity-flow purification. The antibody to be purified should be in 10 to 50 mL of tissue culture supernatant (TCS). Up to 5 mg of antibody can be purified in each run.

This kit is fully compatible with our nanoparticle conjugation kits (Gold, Latex, Europium and Magnetic) (available separately), which allow the purified antibody to be conjugated rapidly.

特記事項 This product is manufactured by Expedeon, an Abcam company. It was previously called AbPure™ TCS Antibody Purification System. Expedeon product code 264-0030 is the same as the 1 pack size of this product and code 264-0500 is the same as the 3 pack size.

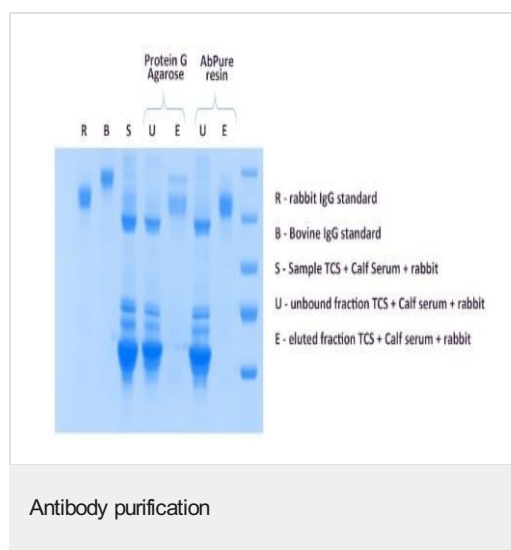
製品の特性

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

内容	1 tests	3 tests
ab273955 - Binding Buffer (10X)	1 x 6ml	1 x 18ml
ab273967 - Collection Tubes (15 mL)	4 units	12 units
ab273953 - Elution Buffer	1 x 7ml	1 x 21ml
ab273956 - Neutralizer	1 x 2ml	1 x 6ml
ab273964 - Preservative	1 x 2ml	1 x 6ml

内容	1 tests	3 tests
Protein A resin	1 x 2ml	3 x 2ml
ab273914 - Spin Cartridge / Collecting Tube Assemblies	1 unit	3 units
ab273966 - Spin Purification Columns	1 unit	3 units
ab273954 - Wash Buffer	1 x 25ml	1 x 75ml

画像



Purification of Rabbit IgG using ab270554.

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