

Product datasheet

Mouse Antibody Purification Kit - Nanoparticles ab204910

Overview

Product name Mouse Antibody Purification Kit - Nanoparticles

Product overview Mouse Antibody Purification Kit - Nanoparticles (ab204910) removes contaminants often found in commercial antibodies (e.g. BSA, glycine, tris and azide) that interfere in conjugation reactions with nanoparticles particles using our nanoparticle conjugation kits (Gold, Latex, Europium and Magnetic). It can also be used to purify mouse antibodies from crude samples such as ascites fluid. The purification system can be used with antibody volumes from 100-500 μ L. 20 to 150 μ g of antibody can be purified in each test.

The method involves capturing the mouse antibody on Mouse Resin and removing unwanted substances by a simple wash procedure, which is performed in a standard microfuge. The purified antibody is then eluted and neutralized.

Antibodies purified using the Mouse Antibody Purification Kit - Nanoparticles are fully compatible with our Gold antibody conjugation kits, Magnetic particle conjugation kits, Latex and Europium conjugation kits (available separately). To purify antibodies for use with our other conjugation kits, we recommend our [Mouse Antibody Purification Kit ab128745](#).

Notes:

(1) The Mouse Antibody Purification Kit - Nanoparticles is not suitable for use with antibodies from other species.

(2) The binding strength of bovine IgG to the Mouse Resin is negligible.

Notes This product is manufactured by Expedeon, an Abcam company, and has previously been called AbPure™ Mouse Specific Purification System and GOLD mouse antibody purification kit. 261-0005 is the same as the 1 test size. 261-0010 is the same as the 3 tests size.

Properties

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

Components	1 tests	3 tests
ab273959 - Additional Collecting Tubes	4 units	12 units
ab273955 - Binding Buffer (10X)	1 x 250µl	1 x 250µl
ab273953 - Elution Buffer	1 x 1500µl	1 x 1500µl
ab273960 - Mouse Resin	1 x 400µl	3 x 400µl
ab273956 - Neutralizer	1 x 250µl	1 x 250µl
ab273958 - Spin cartridge/ collecting tube assemblies	1 unit	3 units
ab273954 - Wash Buffer	1 x 12ml	1 x 12ml

Applications

Our [Abpromise guarantee](#) covers the use of **ab204910** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
AP		Use at an assay dependent concentration.

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