

Product datasheet

Antibody Concentration And Clean-Up Kit – Gold And Magnetic Nanoparticles ab204911

Overview

Product name	Antibody Concentration And Clean-Up Kit – Gold And Magnetic Nanoparticles
Product overview	<p>Antibodies are sometimes only available at concentrations that are too low for antibody conjugation or other antibody applications. The Antibody Concentration And Clean-Up Kit – Gold And Magnetic Nanoparticles uses a simple spin column to easily and quickly remove excess buffer from the antibody thereby providing a more concentrated antibody solution. The kit can also be used to reduce the concentration of additives often found in antibody formulations, such as salt, citrate and azide, that may interfere in subsequent conjugation reactions with our Gold and Magnetic Conjugation Kits. The Antibody Concentration And Clean-Up Kit – Gold And Magnetic Nanoparticles also allows the researcher to perform a simple buffer exchange to transfer the antibody into a more favourable buffer for conjugation using our nanoparticle conjugation kits.</p> <p>Antibodies purified using the Antibody Concentration And Clean-Up Kit – Gold And Magnetic Nanoparticles are fully compatible with our Gold and Magnetic conjugation kits. Please note that kit ab269965 should be used if the antibody is to be used with the Latex and Europium conjugation kits.</p>

Notes	This product is manufactured by Expedeon, an Abcam company, and has previously been called AbPure™ Antibody Concentration and Clean-Up Kit and GOLD antibody concentration kit. 262-0010 is the same as the 3 tests size.
--------------	---

Properties

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

Components	3 tests
ab273914 - Spin Cartridge / Collecting Tube Assemblies	1 unit
ab273961 - Conjugation Buffer	1 x 15ml

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