

Product datasheet

BSA Removal Kit - Nanoparticles ab204912

1 References

Overview

Product name BSA Removal Kit - Nanoparticles

Product overview Our BSA Removal Kit - Nanoparticles (ab204912) is a simple one-step, 10 minute method that effectively separates BSA from the antibody. If necessary, the antibody may then be transferred to a buffer more suited to nanoparticle conjugation. The BSA Removal Kit - Nanoparticles can be used on any antibody sub-type and species. Bovine Serum Albumin (BSA) is often added as a stabilizer to purified antibodies. However, the presence of BSA is not suitable for all antibody applications. For example, during antibody labelling reactions, BSA in the antibody formulation directly competes with the antibody, greatly reducing the conjugation efficiency. It is therefore essential to remove BSA prior to antibody labelling techniques. Common commercial BSA removal techniques can involve many laborious steps.

The BSA Removal Kit can separate BSA from antibody solutions with antibody concentrations from 0.03 mg/mL to 10 mg/mL. Separation is more efficient at higher antibody concentrations. BSA can be effectively separated when present at concentrations of up to 0.5%. If BSA is present at higher concentrations, dilute the antibody mix with de-ionised, distilled water until the BSA concentration is 0.5% or less. 50 µg of antibody is the lower limit for seeing a clearly visible pellet.

Antibodies purified using the BSA Removal Kit - Nanoparticles are fully compatible with our Gold, Magnetic, Latex and Europium conjugation kits (available separately). The BSA Removal Kit can be used on any antibody sub-type, and species. To remove BSA from antibodies prior to conjugation with our other conjugation kits, we recommend using our BSA removal kit ([ab173231](#)).

Important considerations:

The BSA Removal Kit can separate BSA from antibody solutions with antibody concentrations from 0.03 mg/mL to 10 mg/mL. Separation is more efficient at higher antibody concentrations. 50 µg of antibody is the lower limit for seeing a clearly visible pellet.

BSA can be effectively separated when present at concentrations of up to 0.5%. If BSA is present at higher concentrations, dilute the antibody mix with de-ionized, distilled water until BSA concentration is 0.5% or less. Alternatively, if BSA is over a 0.5%, two or more runs may need to be performed to completely remove BSA.

Glycerol concentration must not exceed 20%.

The components of the removal buffer may precipitate at low temperatures. Gently warming the solution should allow solubilisation. Any undissolved crystals should be spun down by brief centrifugation, and the supernatant should be used for the BSA removal.

Notes This product is manufactured by Expedeon, an Abcam company, and has previously been called AbPure™ BSA Removal Kit and GOLD BSA removal kit. 263-0100 is the same as the 1 mL size.

Properties

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

Components	1 ml
ab273962 - BSA Removal Buffer	1 x 1ml
ab273963 - Re-suspension Buffer	1 x 500µl

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