

BSA Removal Kit ab173231

★★★★★ [2 Abreviews](#) [1 图像](#)

概述

产品名称	BSA Removal试剂盒
检测类型	Direct
检测时间	0h 10m
产品概述	<p>Abcam's BSA Removal Kit (ab173231) is a simple one-step, 10 minute method which effectively separates the BSA from the antibody. The antibody is left in a suitable position for transfer to a buffer more suited to conjugation. The BSA Removal Kit can be used on any antibody sub-type, and species.</p>

This 1 ml kit can remove all of the BSA from up to 1.25 ml of antibody, with a BSA concentration of 0.5% or less. For higher BSA concentrations, the method may need to be repeated, or a higher volume of BSA Removal Buffer may be required.

Please note that this kit is not compatible with our Gold, Latex, Europium and Magnetic Conjugation Kits. To remove BSA from antibodies prior to conjugation with these kits, please use BSA Removal Kit - Nanoparticles ([ab204912](#)).

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.

说明

This product is manufactured by Expedeon, an Abcam company, and was previously called AbSelect BSA Removal Kit. 820-0100 is the same as the 1 ml size.

Bovine Serum Albumin (BSA) is often added to purified antibodies as it is an effective stabilizer. However, when labelling antibodies, the BSA becomes a hindrance, as it directly competes with the antibody to attach to the label, greatly reducing the conjugation efficiency. Therefore, prior to undertaking labelling techniques, it is essential to remove the BSA. Common commercial BSA removal techniques can involve many laborious steps.

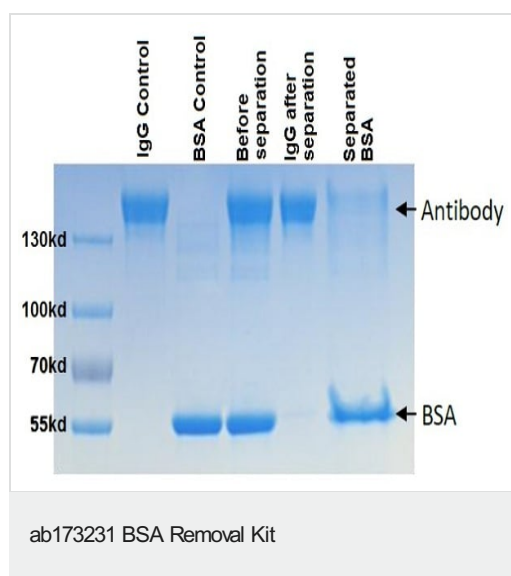
性能

存放说明

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

组件	1 ml
BSA Removal Buffer	1 x 1ml
Re-suspension Buffer	1 x 500µl

图片



SDS-PAGE Gel showing the use of ab173231 BSA Removal Kit on a mixture containing 1 mg/ml IgG and 1 mg/ml BSA. The gel shows the mix before and after separation. 4-12% Bis-Tris gel, non-reducing.

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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors