

# Antibody Purification Kit (Gold/Europium labelling) ab204909

## 概述

### 产品名称

抗体Purification试剂盒(Gold/Europium labelling)

### 产品概述

Antibody Purification Kit - Nanoparticles (ab204909) removes common buffer constituents found in commercial antibodies (e.g. BSA, glycine, tris and azide) that decrease the efficiency of conjugation reactions with nanoparticles using our nanoparticle conjugation kits. ab204909 can also be used to purify antibodies from crude samples such as ascites fluid or serum. The antibody should ideally be in a volume of 100 - 500  $\mu$ L. Up to 500  $\mu$ g of antibody can be purified in each run.

The method involves capture of the antibody on Protein A Resin and removal of unwanted substances by a simple wash procedure, using a standard microfuge. The purified antibody is then eluted and neutralized.

Antibodies purified using the Antibody Purification Kit - Nanoparticles are fully compatible with our nanoparticle conjugation kits (Gold, Latex, Europium and Magnetic). To purify antibodies for use with our other antibody conjugation kits, we recommend using one of our other purification kits - view our [selection guide](#).

For purifying mouse IgG, we recommend using Mouse Antibody Purification Kit - Nanoparticles [ab204910](#).

### 说明

This product is manufactured by Expedeon, an Abcam company, and has previously been called AbPure™ Protein A Purification System and GOLD antibody purification kit. 260-0005 is the same as the 1 test size. 260-0010 is the same as the 3 tests size.

## 性能

### 存放说明

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

组件	1 tests	3 tests
ab273959 - Additional Collecting Tubes	4 vials	12 vials
ab273955 - Binding Buffer (10X)	1 x 250 $\mu$ l	1 x 250 $\mu$ l
ab273953 - Elution Buffer	1 x 1500 $\mu$ l	1 x 1500 $\mu$ l

组件	1 tests	3 tests
ab273956 - Neutralizer	1 x 250µl	1 x 250µl
Protein A resin	1 x 100µl	3 x 100µl
ab273958 - Spin cartridge/ collecting tube assemblies	1 unit	3 units
ab273954 - Wash Buffer	1 x 12ml	1 x 12ml

**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