

Serum Antibody Purification Kit (Protein A) ab109209

1 References

Overview

Product name Serum Antibody Purification Kit (Protein A)

Product overview **Serum Antibody Purification Kit (Protein A) (ab109209)** is prepared by coupling highly purified protein A to agarose beads and can therefore be used to purify IgG fractions from both serum and ascites fluid.

The antibody is captured on the resin and unwanted substances are removed by a simple wash procedure. The purified product is then eluted and neutralized.

The components of **ab109209** are fully compatible with our Lightning-Link® Antibody Conjugation Kits and our Oligonucleotide Conjugation Kit. However they are not compatible with our nanoparticle Antibody conjugation kits (Gold, Latex, Europium and Magnetic). To purify antibodies for use with our nanoparticle conjugation kits, please use our Antibody Purification Kit - Nanoparticles (**ab204909**).

ab109209 is not suitable for goat antibody purification. Please refer to the protocol booklet for more information about suitable and unsuitable species antibodies.

Notes This product is manufactured by Expedeon, an Abcam company, and was previously called AbSelect™ A Serum Antibody Purification System. 863-0500 is the same as the 3 tests size. 863-0030 is the same as the 1 test size.

ab109209 contains 1-3 columns, and each can purify up to **20mg** of antibody per run.

The volume of sample required will depend on the host species (See Table below):

Species	Normal range IgG(mg/ml)	Suitable Volumes for Product(ml)
Rabbit	12 - 15	1.3 - 1.7
Human	7 - 23	0.9 - 2.9
Mouse	2 - 5	4 - 10
Sheep/Goat	18 - 24	0.8 - 1.1
Rat	5 - 7	2.9 - 4

Ascites Fluid	0.5 - 5	4 - 40
---------------	---------	--------

Properties

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

Components	3 tests	1 tests
Binding Buffer (10X)	1 x 18ml	1 x 6ml
Elution Buffer	1 x 15ml	1 x 5ml
Neutralizer Buffer	1 x 2ml	1 x 2ml
<u>ab270308 - Protein A resin</u>	3 x 5ml	1 x 5ml
ab274180 - Purification Column	3 units	1 unit
Wash Buffer	1 x 75ml	1 x 25ml

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