

Product datasheet

Endotoxin Removal Kit (Rapid) ab239707

1 Image

Overview

Product name Endotoxin Removal Kit (Rapid)

Product overview Endotoxin Removal Kit (Rapid)(ab239707) can quickly and effectively eliminate endotoxins to < 0.05 EU/ml in solutions containing proteins or pharmacologically important components via the immobilized polymyxin B, which is known for capturing endotoxin and preventing toxic effects.

The kit quickly and effectively eliminates endotoxins to < 0.05 EU/ml and is ideal for processing small-scale solution sample (0.1-0.5 mL).

Notes This product is manufactured by BioVision, an Abcam company and was previously called K2501 ToxOut™ Rapid Endotoxin Removal Kit. K2501-5 is the same size as the 5 kit size of ab239707.

Properties

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

Components	5 kit
Endotoxin Removal Equilibration Buffer	1 x 10ml
Endotoxin-free Collection Tube	10 units
Rapid Endotoxin Removal Regeneration Buffer	2 x 10ml
Rapid Endotoxin Removal Spin Column	5 units
Rapid Removal Wash Buffer	1 x 10ml

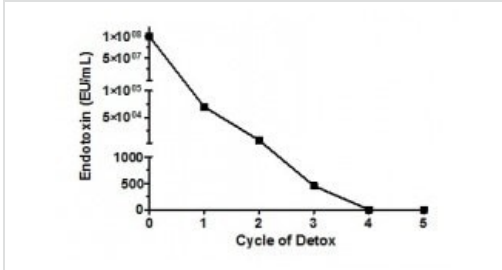
Relevance Endotoxins are part of the outer membrane of the cell wall of gram negative bacteria. Endotoxins are invariably associated with gram negative bacteria whether the organisms are pathogens or not. Although the term "endotoxin" is occasionally used to refer to any cell associated bacterial toxin, it is properly reserved to refer to the lipopolysaccharide complex associated with the outer membrane of gram-negative bacteria such as E. coli, Salmonella, Shigella, Pseudomonas, Neisseria, Haemophilus, and other leading pathogens. The biological activity of endotoxins are associated with lipopolysaccharide (LPS). Toxicity is associated with the lipid component (Lipid A) and immunogenicity is associated with the polysaccharide components. The cell wall antigens (O antigens) of gram negative bacteria are components of LPS. LPS elicits a variety of

inflammatory responses in an animal. Because it activates complement by the alternative (properdin) pathway, it is often part of the pathology of gram negative bacterial infections.

Cellular localization

Cell Membrane

Images



Endotoxin Removal Kit (Rapid)(ab239707) removes >90% of endotoxin from a protein solution sample.

Endotoxin capacities and endotoxin efficiencies were determined by challenging 0.1 mL resin with 1×10^8 EU/mL LPS in 200 μ L BSA (10 mg/mL). By reloading samples to the repeatedly regenerated column, the endotoxin spike is reduced to <0.05 EU/mL in the BSA solution. The highest endotoxin capacity is 9.99×10^8 EU/mL from the first cycle of detox. The average efficiency of 5 cycles of detox is 94.5% and the average protein recovery from 5 cycles of detox is 89.9%.

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