

Datasheet: PUR028

Description:	PROTEUS MIDI ENDOTOXIN REMOVAL STANDARD SPIN COLUMN PACK
Name:	ENDOTOXIN REMOVAL STANDARD KIT
Format:	MIDI - Spin Columns
Product Type:	Purification Kit
Quantity:	12 UNITS

Product Details

Applications

This product has been reported to work in the following applications. This information is derived from testing within our laboratories, peer-reviewed publications or personal communications from the originators. Please refer to references indicated for further information. For general protocol recommendations, please visit www.bio-rad-antibodies.com/protocols.

	Yes	No	Not Determined	Suggested Dilution
Protein Purification	▪			

Where this product has not been tested for use in a particular technique this does not necessarily exclude its use in such procedures.

Preservative Stabilisers

94% Preservative base of resin containing:
20% Aqueous Ethanol
80% 0.1M Sodium Chloride

Endotoxin Level

This kit can be used for samples with endotoxin loads < 30,000 EU. For endotoxin loads < 3,000 EU the Mini Endotoxin Removal kit is recommended. Endotoxin loads up to 1,000,000 EU require the use of a Midi High Capacity Endotoxin Removal kit.

Product Information

This kit offers rapid, one step, high grade (>95%) removal of endotoxin from a wide range of protein solutions, with >90% protein recovery.

We recommend reading the [Proteus Endotoxin Removal Kit Handbook](#) prior to use.

Reagents In The Kit

12 x Spin Column Plugs containing Endotoxin Binding Resin
12 x Midi Standard Spin Columns (20ml capacity)
12 x 50ml Centrifuge Tubes
Plug Insertion Tool
Instructions for use

Additional Reagents Required

0.2µm and 1.2µm Syringe Filter Units
Low Endotoxin Pre-equilibration Buffer e.g PBS pH7.4

Instructions For Use

Instructions for use can be found at www.bio-rad-antibodies.com/uploads/IFU/PUR028.pdf

References

1. Negus, D. *et al.* (2015) Parenteral Administration of Capsule Depolymerase EnvD Prevents Lethal Inhalation Anthrax Infection. [Antimicrob Agents Chemother. pii: AAC.01547-15.](#)

Storage

Upon receipt of kit, remove resin plugs box and store this at +4°C.
DO NOT FREEZE RESIN PLUGS OR STORE AT ROOM TEMPERATURE.

Store remainder of kit at room temperature.

Shelf Life Please see label for expiry date.

Health And Safety Information Material Safety Datasheet documentation available at:
Not required

Regulatory For research purposes only

North & South Tel: +1 800 265 7376

America Fax: +1 919 878 3751

Email: antibody_sales_us@bio-rad.com

Worldwide

Tel: +44 (0)1865 852 700

Fax: +44 (0)1865 852 739

Email: antibody_sales_uk@bio-rad.com

Europe

Tel: +49 (0) 89 8090 95 21

Fax: +49 (0) 89 8090 95 50

Email: antibody_sales_de@bio-rad.com

'M317321:180625'

Printed on 25 Jun 2018

© 2018 Bio-Rad Laboratories Inc | [Legal](#) | [Imprint](#)