

Datasheet: PUR002

Description:	PROTEUS PROTEIN A MIDI PURIFICATION SPIN COLUMN PACK
Name:	ANTIBODY PURIFICATION KIT PROTEIN A
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. Suggested working dilutions are given as a guide only. It is recommended that the user titrates the product for use in their own system using appropriate negative/positive controls.

Preservative Stabilisers

0.01% Thiomersal

Product Information

See [Binding affinities for Protein A and G](#)

Recommended Reading - [Proteus Protein A Antibody Purification Handbook](#)

Reagents In The Kit

12 x Protein A Spin Column Plugs
12 x Spin Columns
24 x 50 ml Centrifuge Tubes
Plug Insertion Tool
Instructions

Storage

Store at +4°C. DO NOT FREEZE.

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

Related Products

Recommended Useful Reagents

[MOUSE MONOCLONAL ANTIBODY ISOTYPING TEST KIT \(MMT1\)](#)

[RAT MONOCLONAL ANTIBODY ISOTYPING TEST KIT \(RMT1\)](#)
[MYCOPLASMA REMOVAL AGENT \(BUF035\)](#)
[PROTEUS PROTEIN G MIDI PURIFICATION SPIN COLUMN PACK \(PUR011\)](#)

North & South America	Tel: +1 800 265 7376 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
----------------------------------	---	------------------	---	---------------	---

'M317328:180625'

Printed on 25 Jun 2018

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