



Datasheet: PUR011

Description:	PROTEUS PROTEIN G MIDI PURIFICATION SPIN COLUMN PACK
Name:	ANTIBODY PURIFICATION KIT PROTEIN G
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

20% Ethanol

Product Information

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

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

Reagents In The Kit

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

References

1. Montaudon, E. *et al.* (2015) Cardiac effects of long-term active immunization with the second extracellular loop of human β 1- and/or β 3-adrenoceptors in Lewis rats. [Pharmacol Res. 100: 210-9](#).

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 A MIDI PURIFICATION SPIN COLUMN PACK \(PUR002\)](#)

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

'M317323:180625'

Printed on 25 Jun 2018

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