

Datasheet: PUR007

Description:	PROTEUS PROTEIN A MINI PURIFICATION SPIN COLUMN PACK
Name:	ANTIBODY PURIFICATION KIT PROTEIN A
Format:	MINI - Spin Columns
Product Type:	Purification Kit
Quantity:	48 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

48 x Protein A Spin Column Plugs
48 x Spin Columns
96 x 2.2 ml Centrifuge Tubes
Plug Insertion Tool
Instructions

References

1. Bekesova, S. *et al.* (2012) N-glycans in liver-secreted and immunoglobulin-derived protein fractions. [J Proteomics. 75 \(7\): 2216-24.](#)
2. Goldman, R. (2012) Glycomic Analysis of Prostate Cancer [Medicine and Medical Research Award Number: W81XWH-09-1-0202](#)

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 MINI PURIFICATION SPIN COLUMN PACK \(PUR015\)](#)

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

'M317329:180625'

Printed on 25 Jun 2018

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