

Datasheet: PUR012

Description:	PROTEUS PROTEIN G MIDI PURIFICATION KIT
Name:	ANTIBODY PURIFICATION KIT PROTEIN G
Format:	MIDI - Spin Columns, Buffers, U/F Spinners
Product Type:	Purification Kit
Quantity:	4 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

Buffers contain 0.09% Sodium Azide (NaN₃)
Resin contains 20% Ethanol

Product Information

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

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

Reagents In The Kit

4 x Protein G Spin Column Plugs
4 x Spin Columns
8 x 50 ml Centrifuge Tubes
4 x 10 kDa MWCO Ultrafiltration Spinners
1 x 250 ml of 1 x Buffer G
1 x 125 ml of 1 x Buffer B2
1 x 30 ml of 1 x Buffer C
Plug Insertion Tool
Instructions

References

1. Abdelkrim, M.A. *et al.* (2012) Autoantibodies against cardiac $\beta(1)$ -adrenoceptor do not affect the low-affinity state $\beta(1)$ -adrenoceptor-mediated inotropy in rat cardiomyocytes. [Can J Physiol Pharmacol. 90: 407-14.](#)
2. Soler, L. *et al.* (2011) Development and validation of a novel competitive ELISA for the detection of serum amyloid A in pigs. [Vet J. 190: e7-11.](#)
3. Kimura, K. *et al.* (2012) Pyrroloquinoline quinone stimulates epithelial cell proliferation by activating epidermal growth factor receptor through redox cycling. [Free Radic Biol Med. 53: 1239-51.](#)
4. Montaudon, E. *et al.* (2014) Cardiovascular effects of beta1 and beta3-adrenergic receptor autoantibodies in Lewis rat.

[Article in French] [Ann Cardiol Angeiol \(Paris\). 63: 128-34.](#)

5. Yuan, W. *et al.* (2015) Quantitative analysis of immunoglobulin subclasses and subclass specific glycosylation by LC-MS-MRM in liver disease. [J Proteomics. 116: 24-33.](#)

6. Broecker, F. *et al.* (2016) Multivalent display of minimal *Clostridium difficile* glycan epitopes mimics antigenic properties of larger glycans. [Nat Commun. 7: 11224.](#)

7. 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 #10125 #10253 #10254 available at: Protein G Binding Buffer G (10125): https://www.bio-rad-antibodies.com/uploads/MSDS/10125.pdf Elution Buffer B2 (10253): https://www.bio-rad-antibodies.com/uploads/MSDS/10253.pdf Neutralisation Buffer C (10254): https://www.bio-rad-antibodies.com/uploads/MSDS/10254.pdf
--------------------------------------	---

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 KIT \(PUR003\)](#)

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

'M317320:180625'

Printed on 25 Jun 2018

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