

Datasheet: PUR016

Description:	PROTEUS PROTEIN G MINI PURIFICATION KIT
Name:	ANTIBODY PURIFICATION KIT PROTEIN G
Format:	MINI - Spin Columns, Buffers, U/F Spinners
Product Type:	Purification Kit
Quantity:	16 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 (NaN ₃) 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	16 x Protein G Spin Column Plugs 16 x Spin Columns 32 x 2.2 ml Centrifuge Tubes 16 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

- 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.](#)
- 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.](#)
- 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.](#)
- 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 MINI PURIFICATION KIT \(PUR008\)](#)

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

'M317325:180625'

Printed on 25 Jun 2018

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