

Datasheet: PUR008

Description:	PROTEUS PROTEIN A MINI PURIFICATION KIT
Name:	ANTIBODY PURIFICATION KIT PROTEIN A
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 (NaN₃)
Resin contains 0.01% Thiomersal

Product Information

See [Binding affinities for Protein A and G](#). Recommended Reading - [Proteus Protein A Purification Handbook](#).

Reagents In The Kit

16 x Protein A 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 A
1 x 125 ml of 1 x Buffer B1
1 x 125 ml of 1 x Buffer B2
1 x 30 ml of 1 x Buffer C
Plug Insertion Tool
Instructions

References

1. Golchin, M. *et al.* (2012) Latex agglutination test based on single-chain Fv recombinant antibody fragment. [Scand J Immunol. 75: 38-45.](#)
2. 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.](#)
3. Jansen, C.A. *et al.* (2016) Identification of an Activating Chicken Ig-like Receptor Recognizing Avian Influenza Viruses. [J Immunol. 197 \(12\): 4696-703.](#)

Storage

Store at +4°C. DO NOT FREEZE.

Shelf Life Please see label for expiry date.

Health And Safety Information Material Safety Datasheet documentation #10251 #10253 #10254 #10252 available at:
Protein A Binding Buffer A (10251): <https://www.bio-rad-antibodies.com/uploads/MSDS/10251.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>
Elution Buffer B1 (10252): <https://www.bio-rad-antibodies.com/uploads/MSDS/10252.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 G MINI PURIFICATION KIT \(PUR016\)](#)

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
----------------------------------	---	------------------	---	---------------	---

'M317322:180625'

Printed on 25 Jun 2018

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