

Datasheet: PUR014

Description:	PROTEUS PROTEIN G MINI PURIFICATION STARTER KIT
Name:	ANTIBODY PURIFICATION KIT PROTEIN G
Format:	MINI - Spin Columns, Buffers, U/F Spinners
Product Type:	Purification Kit
Quantity:	2 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	2 x Protein G Spin Column Plugs 2 x Spin Columns 4 x 2.2 ml Centrifuge Tubes 2 x 10 kDa MWCO Ultrafiltration Spinners 1 x 60 ml of 1 x Buffer G 1 x 20 ml of 1 x Buffer B2 1 x 10 ml of 1 x Buffer C Plug Insertion Tool Instructions
----------------------------	---

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 \(RMT1\)](#)

[PROTEUS PROTEIN A MINI PURIFICATION STARTER KIT \(PUR006\)](#)

[MYCOPLASMA REMOVAL AGENT \(BUF035\)](#)

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

'M317324:180625'

Printed on 25 Jun 2018

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