

Datasheet: PUR006

Description:	PROTEUS PROTEIN A MINI PURIFICATION STARTER KIT
Name:	ANTIBODY PURIFICATION KIT PROTEIN A
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 0.01% Thiomersal

Product Information

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

Reagents In The Kit

2 x Protein A 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 A
1 x 20 ml of 1 x Buffer B1
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 #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>

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 STARTER KIT \(PUR014\)](#)

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

'M317319:180625'

Printed on 25 Jun 2018