

Datasheet: PUR023

Description:	PROTEUS IMAC MINI PURIFICATION SPIN COLUMN PACK + BUFFERS
Name:	HIS-TAG PROTEIN
Format:	MINI - Spin Columns, Buffers
Product Type:	Purification Kit
Quantity:	24 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	IMAC Ni-IDA binds to HIS-6 tagged proteins.
----------------------------	---

Recommended Reading - [Proteus IMAC Purification Handbook](#).

Reagents In The Kit	24 x Ni-IDA Spin Column Plugs 24 x Spin Columns 48 x 2.2 ml Centrifuge Tubes 1 x 250 ml of 5 x Buffer A 1 x 125 ml of 1 x Buffer B Plug Insertion Tool Instructions
----------------------------	---

Storage	Store at +4°C. DO NOT FREEZE.
----------------	-------------------------------

Shelf Life	24 months from date of despatch.
-------------------	----------------------------------

Health And Safety Information	Material Safety Datasheet documentation available at: Material Safety Datasheet Documentation #10264, #10156 available at: https://www.bio-rad-antibodies.com/uploads/MSDS/10264.pdf https://www.bio-rad-antibodies.com/uploads/MSDS/10156.pdf
--------------------------------------	---

Regulatory	For research purposes only
-------------------	----------------------------

Related Products

Recommended Useful Reagents

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

[MOUSE ANTI HISTIDINE TAG \(MCA1396\)](#)

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

'M310359:170919'

Printed on 19 Sep 2017

© 2017 Bio-Rad Laboratories Inc | [Legal](#) | [Imprint](#)