

Datasheet: PUR019

Description:	PROTEUS IMAC MIDI PURIFICATION SPIN COLUMN PACK + BUFFERS
Name:	HIS-TAG PROTEIN
Format:	MIDI - Spin Columns, Buffers
Product Type:	Purification Kit
Quantity:	8 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 20% Ethanol

Product Information

IMAC Ni-IDA binds to HIS-6 tagged proteins.

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

Reagents In The Kit

8 x Ni-IDA Spin Column Plugs
8 x Spin Columns
12 x 50 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

References

1. Inobe, T. *et al.* (2015) Artificial regulation of p53 function by modulating its assembly. [Biochem Biophys Res Commun. 467 \(2\): 322-7.](#)

Storage

Store at +4°C. DO NOT FREEZE.

Shelf Life

24 months from date of despatch.

Acknowledgements

His-tag is a registered trademark of EMD Biosciences.

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

'M313881:180403'

Printed on 03 Apr 2018

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