

## Datasheet: PUR026

<b>Description:</b>	PROTEUS IMAC MINI PURIFICATION KIT
<b>Name:</b>	HIS-TAG PROTEIN
<b>Format:</b>	MINI - Spin Columns, Buffers, U/F Spinners
<b>Product Type:</b>	Purification Kit
<b>Quantity:</b>	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](http://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.

<b>Preservative Stabilisers</b>	Buffers contain 0.09% Sodium Azide (NaN <sub>3</sub> ) Resin contains 20% Ethanol
---------------------------------	--

<b>Product Information</b>	IMAC Ni-IDA binds to HIS-6 tagged proteins.
----------------------------	---

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

<b>Reagents In The Kit</b>	24 x Ni-IDA Spin Column Plugs 24 x Spin Columns 48 x 2.2 ml Centrifuge Tubes 24 x 10 kDa MWCO Ultrafiltration Spinners 1 x 250 ml of 5 x Buffer A 1 x 125 ml of 1 x Buffer B Plug Insertion Tool Instructions
----------------------------	--

<b>References</b>	<ol style="list-style-type: none"> <li>Lombardi, A. <i>et al.</i> (2010) Pichia pastoris as a host for secretion of toxic saporin chimeras. <a href="#">FASEB J. 24: 253-65.</a></li> <li>Inobe, T. <i>et al.</i> (2015) Artificial regulation of p53 function by modulating its assembly. <a href="#">Biochem Biophys Res Commun. 467 (2): 322-7.</a></li> <li>Inobe, T. &amp; Nozaki, M. (2016) Proteasomal degradation of damaged polyubiquitin. <a href="#">Biochem Biophys Res Commun. pii: S0006-291X(16)30196-6. [Epub ahead of print]</a></li> <li>Inobe, T. &amp; Nukina, N. (2016) Rapamycin-induced oligomer formation system of FRB-FKBP fusion proteins. <a href="#">J Biosci Bioeng. 122 (1): 40-6.</a></li> </ol>
-------------------	--

<b>Further Reading</b>	1. <a href="#">Proteus IMAC Purification Handbook</a> . (2012) Purification Handbook
------------------------	--

<b>Storage</b>	Store at +4°C. DO NOT FREEZE.
<b>Shelf Life</b>	24 months from date of despatch.
<b>Health And Safety Information</b>	Material Safety Datasheet documentation available at: Material Safety Datasheet Documentation #10264, #10156 available at: <a href="https://www.bio-rad-antibodies.com/uploads/MSDS/10264.pdf">https://www.bio-rad-antibodies.com/uploads/MSDS/10264.pdf</a> <a href="https://www.bio-rad-antibodies.com/uploads/MSDS/10156.pdf">https://www.bio-rad-antibodies.com/uploads/MSDS/10156.pdf</a>
<b>Regulatory</b>	For research purposes only

## Related Products

### Recommended Useful Reagents

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

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

<b>North &amp; South America</b>	Tel: +1 800 265 7376 Fax: +1 919 878 3751 Email: <a href="mailto:antibody_sales_us@bio-rad.com">antibody_sales_us@bio-rad.com</a>	<b>Worldwide</b>	Tel: +44 (0)1865 852 700 Fax: +44 (0)1865 852 739 Email: <a href="mailto:antibody_sales_uk@bio-rad.com">antibody_sales_uk@bio-rad.com</a>	<b>Europe</b>	Tel: +49 (0) 89 8090 95 21 Fax: +49 (0) 89 8090 95 50 Email: <a href="mailto:antibody_sales_de@bio-rad.com">antibody_sales_de@bio-rad.com</a>
----------------------------------	---	------------------	---	---------------	---

'M310361:170919'

**Printed on 19 Sep 2017**