

## Datasheet: PUR025

<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>	4 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.

### Preservative Stabilisers

Buffers contain 0.09% Sodium Azide (NaN<sub>3</sub>)  
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

4 x Ni-IDA Spin Column Plugs  
4 x Spin Columns  
8 x 2.2 ml Centrifuge Tubes  
4 x 10 kDa MWCO Ultrafiltration Spinners  
1 x 60 ml of 5 x Buffer A  
1 x 35 ml of 1 x Buffer B  
Plug Insertion Tool  
Instructions

### References

- Lombardi, A. *et al.* (2010) Pichia pastoris as a host for secretion of toxic saporin chimeras. [FASEB J. 24: 253-65.](#)
- Inobe, T. *et al.* (2015) Artificial regulation of p53 function by modulating its assembly. [Biochem Biophys Res Commun. 467 \(2\): 322-7.](#)
- Inobe, T. & Nozaki, M. (2016) Proteasomal degradation of damaged polyubiquitin. [Biochem Biophys Res Commun. pii: S0006-291X\(16\)30196-6. \[Epub ahead of print\]](#)
- Inobe, T. & Nukina, N. (2016) Rapamycin-induced oligomer formation system of FRB-FKBP fusion proteins. [J Biosci Bioeng. 122 \(1\): 40-6.](#)

### 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 #10264 #10156 available at:  
PBS Buffer A (10264): <https://www.bio-rad-antibodies.com/uploads/MSDS/10264.pdf>  
Imidazole Buffer B (10156): <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](mailto: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](mailto: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](mailto:antibody_sales_de@bio-rad.com)

'M317331:180625'

**Printed on 25 Jun 2018**

---

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