

Datasheet: BUF039A

Description:	ELISA NEPTUNE ASSAY DILUENT
Name:	ELISA NEPTUNE ASSAY DILUENT
Format:	Ready To Use
Product Type:	Accessory Reagent
Quantity:	100 ml

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
ELISA	▪			Neat

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.

Product Form	Ready to use - liquid
Buffer Solution	Phosphate buffered saline
Preservative Stabilisers	0.1% Sodium Azide
Instructions For Use	Pipette buffer directly onto the plate just prior to adding the sample or standard (suggested volume 50 - 100ul/well).
References	<ol style="list-style-type: none"> Defresne, F. <i>et al.</i> (2010) Differential influence of anticancer treatments and angiogenesis on the seric titer of autoantibody used as tumor and metastasis biomarker. Neoplasia 12: 562-570. Massei, G. <i>et al.</i> (2015) Immunocontraception for managing feral cattle in Hong Kong. PLoS One. 10 (4): e0121598.
Storage	Store at +4°C. DO NOT FREEZE
Shelf Life	18 months from date of despatch.
Health And Safety Information	Material Safety Datasheet documentation #10377 available at: 10377: https://www.bio-rad-antibodies.com/uploads/MSDS/10377.pdf
Regulatory	For research purposes only

Related Products

Recommended Useful Reagents

[ELISA ULTRABLOCK \(BUF033A\)](#)

[ELISA ULTRABLOCK \(BUF033B\)](#)

[ELISA ULTRABLOCK \(BUF033C\)](#)

[5x ELISA COATING BUFFER \(BUF030A\)](#)

[10x ELISA WASH BUFFER \(BUF031A\)](#)

[ELISA BSA BLOCK \(BUF032A\)](#)

North & South America Tel: +1 800 265 7376
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

'M308736:170810'

Printed on 30 Apr 2018

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