

Datasheet: BUF038A

Description:	ELISA IgM REDUCING ASSAY DILUENT
Name:	ELISA IgM REDUCING 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	0.10% Sodium Azide
Stabilisers	1-5% Sodium Metabisulphite

Intended Use

BUF038A is an ELISA assay diluent designed to reduce background noise and assist in matrix equalization in problematic serum and plasma samples, especially those affected by high levels of “sticky” IgM which can non-specifically bind the HRP conjugate. Specially formulated for use in sandwich ELISA, it contains mammalian protein additives based on normal goat serum to reduce non-specific interactions between sample proteins and the plate surface.

The difference in complexity, between serum or plasma samples and diluted standard curve proteins, is a potential cause of target analyte under-recovery. BUF038A equalizes the matrix, reduces IgM-mediated bridging interference, inhibits complement and thrombin activity and contains an antimicrobial agent.

Instructions For Use	Pipette buffer directly onto the plate, just prior to adding the sample or standard (suggested volume 50 - 100ul/well).
	NB. Since this buffer contains proteins derived from goat, BUF038A is not recommended for ELISA where an anti goat secondary antibody is used.

Storage	Store at +4°C.
	DO NOT FREEZE

Shelf Life	Please see label for expiry date.
Health And Safety Information	Material Safety Datasheet documentation available at: Material Safety Datasheet Documentation #10376 available at: https://www.bio-rad-antibodies.com/uploads/MSDS/10376.pdf
Regulatory	For research purposes only

Related Products

Recommended Useful Reagents

[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
----------------------------------	---	------------------	---	---------------	---

'M312917:180305'

Printed on 05 Mar 2018

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