



Datasheet: BUF038B

Description:	ELISA IgM REDUCING ASSAY DILUENT
Name:	ELISA IgM REDUCING ASSAY DILUENT
Format:	Ready To Use
Product Type:	Accessory Reagent
Quantity:	500 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 BUF038B 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. BUF038B 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, BUF038B is not recommended for ELISA where an anti goat secondary antibody is used.

Storage Store at +4°C.

DO NOT FREEZE

Shelf Life	18 months from date of despatch.
-------------------	----------------------------------

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 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

'M262379:141216'

Printed on 05 Mar 2018

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