

Datasheet: BUF037C

Description:	ELISA GENERAL ASSAY DILUENT
Name:	ELISA GENERAL ASSAY DILUENT
Format:	Ready To Use
Product Type:	Accessory Reagent
Quantity:	1000 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

0.17% Proclin™ 300

Intended Use

BUF037C is an ELISA assay diluent designed to reduce background noise and assist in matrix equalization in problematic serum, plasma or cell culture samples. Designed for use in all sandwich ELISA formats, it contains mammalian serum proteins 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. BUF037C equalizes the matrix, 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 and standard (suggested volume 50 - 100 ul/well).

NB. Since this buffer contains proteins derived from goat, BUF037C 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.

Acknowledgements Proclin™ 300 is a trademark of The Dow Chemical Company (“Dow”) or an affiliated company of Dow.

Health And Safety Information Material Safety Datasheet documentation #10375 available at:
10375: <https://www.bio-rad-antibodies.com/uploads/MSDS/10375.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

'M295044:161013'

Printed on 10 May 2018

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