



Datasheet: OBT1998GX

Description:	ELISA PLASMA SAMPLE DILUENT
Name:	ELISA PLASMA SAMPLE 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
<10% Bovine Serum Albumin

Intended Use

OBT1998GX is a sample diluent specially formulated for use with serum and plasma samples in antigen-down and sandwich ELISA formats. OBT1998GX dilutes highly concentrated samples into the functional range of the assay, preventing overloading of the coating proteins. It also performs well as a protein diluent in standard curve preparation. The proprietary additives in a buffered BSA protein base minimize non-specific binding and inhibit clotting and interference caused by complement activity. An antimicrobial agent provides a solution free of bacterial contamination.

Instructions For Use

Dilute samples as required before addition to ELISA plate. Highly concentrated samples may need to be diluted 1:1000 or more.

Storage

Store at +4°C. DO NOT FREEZE.
This product should be stored undiluted.

Shelf Life

Please see label for expiry date

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 available at:
Material Safety Datasheet Documentation #10388 available at:

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

'M308624:170727'

Printed on 18 Jan 2018