

Datasheet: BUF033C

Description:	ELISA ULTRABLOCK
Name:	ELISA ULTRABLOCK
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 antibody 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 antibody for use in their own system using the appropriate negative/positive controls.

Product Form Ready to use - liquid

Buffer Solution Phosphate buffered saline

Preservative Stabilisers 0.05% Sodium Azide

Intended Use BUF033C is a high performance ELISA blocking buffer for use in all ELISA formats with high background problems. The formulation includes a proprietary fish extract, molecular stabilizers and an antimicrobial agent. It is particularly useful in antigen-down and sandwich ELISA assays with mammalian samples, especially human, bovine and porcine serum, where the non-mammalian blocking proteins are less likely to interact with endogenous antibodies in the serum sample.

Superior blocking is possible due to the small size of the molecules which block non-specific binding sites on the adsorbed protein and unoccupied regions of the polystyrene plates often sterically inaccessible to traditional blockers. It provides a long-term stable environment for coating antigen or capture antibody. Plates can be blocked at room temperature and stored once dried for up to a year at +4°C.

- Instructions For Use**
1. Coat ELISA plate with antibody or antigen as required.
 2. After incubation, remove the coating solution and wash the plate x2 with wash buffer. [BUF031A](#) can be used for this purpose.
 3. Add 300-400ul of BUF033C and incubate for 2-24 hours. Use a volume equal to or greater than the volume of coating solution.

4. After removal of the blocking buffer continue with the assay or dry the plate for long-term storage at +4°C.

References

1. Colwell, D.D. *et al.* (2010) *Dicrocoelium dendriticum* in cattle from Cypress Hills, Canada: Humoral response and preliminary evaluation of an ELISA. [Vet Parasitol.174: 162-165.](#)
2. Defresne, F. *et al.* (2010) Differential influence of anticancer treatments and angiogenesis on the seric titer of autoantibody used as tumor and metastasis biomarker. [Neoplasia 12: 562-570.](#)
3. Zhang, Y. *et al.* (2012) Causes of alternative pathway dysregulation in dense deposit disease. [Clin J Am Soc Nephrol. 7 \(2\): 265-74.](#)
4. Krachudel, J. *et al.* (2013) Luteal insufficiency in bitches as a consequence of an autoimmune response against progesterone? [Theriogenology. 79: 1278-83.](#)

Storage

Store at +4°C.

DO NOT FREEZE

This product should be stored undiluted.

Shelf Life

18 months from date of despatch.

Health And Safety Information

Material Safety Datasheet documentation available at:
Material Safety Datasheet Documentation #10379 available at:
<https://www.bio-rad-antibodies.com/uploads/MSDS/10379.pdf>

Regulatory

For research purposes only

Related Products

Recommended Useful Reagents

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

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

[ELISA NEPTUNE ASSAY DILUENT \(BUF039A\)](#)

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

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

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

'M296234:161102'

Printed on 13 Sep 2017

© 2017 Bio-Rad Laboratories Inc | [Legal](#) | [Imprint](#)