

Datasheet: BUF032A

Description:	ELISA BSA BLOCK
Name:	ELISA BSA BLOCK
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 systems 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

BUF032A is a general purpose ELISA blocking buffer for use in all ELISA formats requiring low to average blocking strength. The buffer contains a mammalian protein blocking agent based on BSA, proprietary protein stabilizers and an antimicrobial agent. It provides a long-term stable environment for coating antigen or capture antibody, whilst minimizing non-specific binding interactions in the assay process. 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 BUF032A 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.

Storage

Store at +4°C.

DO NOT FREEZE

This product should be stored undiluted.

Shelf Life Please see label for expiry date.

Health And Safety Information Material Safety Datasheet documentation available at:
Material Safety Datasheet Documentation #10378 available at:
<https://www.bio-rad-antibodies.com/uploads/MSDS/10378.pdf>

Regulatory For research purposes only

Related Products

Recommended Useful Reagents

[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

'M309016:170817'

Printed on 22 Nov 2017

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