

Datasheet: BUF032C

Description:	ELISA BSA BLOCK
Name:	ELISA BSA BLOCK
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	0.05% Sodium Azide
Intended Use	BUF032C 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 BUF032C 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

'M311429:171122'

Printed on 22 Nov 2017

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