

Datasheet: BUF030B

Description:	5x ELISA COATING BUFFER
Name:	5x ELISA COATING BUFFER
Format:	5 X Concentrate
Product Type:	Accessory Reagent
Quantity:	500 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	■			

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.

Product Form	5 x concentrate - liquid
Buffer Solution	Phosphate buffered saline
Preservative Stabilisers	0.5% Sodium Azide
Intended Use	BUF030B is a specialist coating buffer for use in all ELISA formats. Designed to stabilize the tertiary structure of the antibody or antigen and maximize adsorption onto polystyrene plates. It is especially effective in sandwich ELISA, preserving the antigen recognition regions of the coating antibody, allowing greater binding reactivity. BUF030B contains an antimicrobial agent allowing coating to be performed at room temperature. Dried plates can be kept at 4°C for long term storage.
Instructions For Use	1. Dilute the required volume of BUF030B 1:5 with deionised water and mix for 5-15 minutes to prepare the working solution. 2. Add antibody or antigen to the working solution and mix for 5-15 minutes. 3. Coat plate (suggested volume 100-200µl per well) and incubate at room temperature for 6-24 hours, protected from light and covered to avoid evaporation.
Storage	Store at +4°C. DO NOT FREEZE. This product should be stored undiluted. Should this product contain a precipitate, heat gently until dissolved. Do not boil.
Shelf Life	Please see label for expiry date.

Health And Safety Information Material Safety Datasheet documentation available at:
Material Safety Datasheet Documentation #10383 available at:
<https://www.bio-rad-antibodies.com/uploads/MSDS/10383.pdf>

Regulatory For research purposes only

Related Products

Recommended Useful Reagents

[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

'M312922:180305'

Printed on 05 Mar 2018

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