

## Datasheet: BUF030A

<b>Description:</b>	5x ELISA COATING BUFFER
<b>Name:</b>	5x ELISA COATING BUFFER
<b>Format:</b>	5 X Concentrate
<b>Product Type:</b>	Accessory Reagent
<b>Quantity:</b>	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](http://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.

<b>Product Form</b>	5 x concentrate - liquid
<b>Buffer Solution</b>	Phosphate buffered saline
<b>Preservative Stabilisers</b>	0.5% Sodium Azide

### Intended Use

BUF030A 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.

BUF030A 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 BUF030A 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](mailto: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](mailto: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](mailto:antibody_sales_de@bio-rad.com)

'M310651:180108'

**Printed on 05 Mar 2018**

---

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