

## ELISA Assay Diluent (5X)

<b>Catalog# / Size</b>	421203 / 100 mL
<b>Regulatory Status</b>	RUO
<b>Other Names</b>	ELISA assay diluent, ELISA diluent
<b>Description</b>	ELISA Assay Diluent is a 5x concentrated phosphate buffered saline solution containing (pre-screened) bovine serum. This buffer is specifically formulated and tested for ELISA applications, such as blocking antibody-coated ELISA plate microwells, diluting ELISA standards and samples, and diluting biotinylated detection antibodies or tertiary reagents (specifically Avidin-HRP or Streptavidin-HRP). The diluent contains a preservative compatible with peroxidase mediated enzyme activity.

### Product Details

---

**Storage & Handling** This solution can be stored unopened between 2°C and 8°C. After opening, it is recommended to use within 30 days or to be aliquoted and stored at -20°C (frost-free) or -70°C for long term storage.

Warm ELISA reagents to room temperature (18-25°C) before use.

**Application** [ELISA](#)

**Recommended Usage** Prior to use, dilute the ELISA Assay Diluent (5X) in PBS, pH 7.4.

PBS recipe: 8.0 g NaCl, 1.16 g Na<sub>2</sub>HPO<sub>4</sub>, 0.2 g KH<sub>2</sub>PO<sub>4</sub>, and 0.2 g KCl. Add deionized water to 1.0 L, filter through 0.2 µm filter, pH to 7.4

**Application References**

(PubMed link indicates BioLegend citation)

1. Kianitsa K, *et al.* 2014. *Nucleic Acids Res.* 42:108. [PubMed](#)

**Product Citations**

1. Maizels K 2014. *Nucleic Acids Res.* 42:108. [PubMed](#)

### Antigen Details

---

**Gene ID** NA

### Related Protocols

---

[Sandwich ELISA Protocol](#)

For research use only. Not for diagnostic use. Not for resale. BioLegend will not be held responsible for patent infringement or other violations that may occur with the use of our products.

\*These products may be covered by one or more Limited Use Label Licenses (see the BioLegend Catalog or our website, [www.biolegend.com/ordering#license](http://www.biolegend.com/ordering#license)). BioLegend products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products, reverse engineer functionally similar materials, or to provide a service to third parties without written approval of BioLegend. By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.

BioLegend Inc., 8999 BioLegend Way, San Diego, CA 92121 [www.biolegend.com](http://www.biolegend.com)  
Toll-Free Phone: 1-877-Bio-Legend (246-5343) Phone: (858) 768-5800 Fax: (877) 455-9587