

ELISA Wash Buffer (20X)

Catalog# / Size 421601 / 500 mL
Regulatory Status RUO

Product Details

Formulation 20X concentrated Phosphate buffered saline (pH7.2) with Tween-20.

Storage & Handling The undiluted solution should be stored at room temperature to avoid salt crystallization.

Application [ELISA](#)

Recommended Usage The BioLegend ELISA Wash Buffer (20X) must be diluted to 1X working solution with D.I. water prior to ELISA wash procedures. e.g. Dilute one (1) part ELISA Wash Buffer (20X) to nineteen (19) parts D.I. water. The 1X solution contains a final concentration of 0.05% Tween-20. Each lot of BioLegend ELISA Wash Buffer (20X) has been tested by ELISA.

Application References
(PubMed link indicates BioLegend citation)

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

Product Citations

1. Sinani G, *et al.* 2017. *Int J Pharm.* 10.1016/j.ijpharm.2017.06.053. [PubMed](#)
2. Lim J, *et al.* 2020. *Sci Rep.* 10:14897. [PubMed](#)
3. 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). 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.

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