

ELISA Coating Buffer (5X)

Catalog# / Size	421701 / 30 mL
Regulatory Status	RUO

Product Details

Formulation	5X concentrated carbonate buffer, pH 9.5.
Storage & Handling	The undiluted solution should be stored between 2°C and 8°C.
Application	ELISA - Quality tested
Recommended Usage	The BioLegend ELISA Coating Buffer(5X) must be diluted to 1X working solution with D.I. water prior to ELISA coating procedures. e.g. Dilute one (1) part ELISA Coating Buffer(5X) to four (4) parts D.I. water. Each lot of BioLegend ELISA Coating Buffer (5X) has been tested with cytokine ELISA capture antibodies.

Application References

(PubMed link indicates BioLegend citation)

1. Kianitsa K, *et al.* 2014. *Nucleic Acids Res.* 42:108. [PubMed](#)
2. Fritz JM, *et al.* 2014. *PLoS One.* 9:107473. [PubMed](#)

Product Citations

1. Maizels K 2014. *Nucleic Acids Res.* 42:108. [PubMed](#)
2. Fritz J, Weaver T 2014. *PLoS One.* 9:107473. [PubMed](#)
3. Lakshmi Priya T, *et al.* 2016. *PLoS One.* 11: 0151153. [PubMed](#)
4. Lakshmi Priya T, *et al.* 2017. *Int J Biol Macromol.* 10.1016/j.ijbiomac.2017.07.115. [PubMed](#)
5. Yoon I, *et al.* 2017. *Vaccine.* 10.1016/j.vaccine.2017.07.037. [PubMed](#)
6. Lim J, *et al.* 2020. *Sci Rep.* 10:14897. [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.

BioLegend Inc., 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