



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

Immuno Buffer 1

Catalog #: 79311

Description: The 3X Immuno Buffer 1 is optimized for use with a large variety of Immunotherapy Assay Kits.

Applications: Suitable for studying enzyme kinetics, screening small molecular inhibitors, and for high throughput applications in drug discovery.

Supplied As: Phosphate-buffered saline solution

Format: Aqueous buffer solution

Quantity: 50 ml

Storage/Stability: Store at -20°C. Avoid freeze/thaw cycles. At least 12 months from date of receipt, when stored as directed.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: [**info@bpsbioscience.com**](mailto:info@bpsbioscience.com)
Please visit our website at: [**www.bpsbioscience.com**](http://www.bpsbioscience.com)



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Optimized For Use With:

Product	Cat. #	Size
CD28:B7-1[Biotinylated] Inhibitor Screening Assay Kit	72007	96 rxns
CD28:B7-2[Biotinylated] Inhibitor Screening Assay Kit	72062	96 rxns
CD47:SIRP- α [Biotinylated] Inhibitor Screening Assay Kit	72044	96 rxns
CD47:SIRP- γ [Biotinylated] Inhibitor Screening Assay Kit	72059	96 rxns
CD137[Biotinylated]:CD137L Inhibitor Screening Assay Kit	72025	96 rxns
CD226:CD155 Homogeneous Assay Kit	72052	384 rxns
CD226:CD112 Homogeneous Assay Kit	72051	384 rxns
CTLA4:B7-1 TR-FRET Assay Kit	72120	384 rxns
CTLA4:B7-1[Biotinylated] Inhibitor Screening Assay Kit	72009	96 rxns
CTLA4[Biotinylated]:B7-2 Inhibitor Screening Assay Kit	72024	96 rxns
IL-17RA[Biotin]:IL-17A Inhibitor Screening Assay Kit	72060	96 rxns
OX40[Biotinylated]:OX40L Inhibitor Screening Assay Kit	72045	96 rxns
PD-1:PD-L2[Biotinylated] Inhibitor Screening Colorimetric Assay Kit	72017	96 rxns
PD-1[Biotinylated]:PD-L1 Homogeneous Assay Kit	72028	384 rxns
PD-1[Biotinylated]:PD-L1 Inhibitor Screening Assay Kit	72005	96 rxns
PD-1[Biotinylated]:PD-L1 Inhibitor Screening Colorimetric Assay Kit	72018	96 rxns
PD-1[Biotinylated]:PD-L2 Inhibitor Screening Assay Kit	72006	96 rxns
PD-1[Biotinylated]:PD-L2 Inhibitor Screening Colorimetric Assay Kit	72019	96 rxns
PD-L1:B7-1[Biotinylated] Inhibitor Screening Assay Kit	72026	96 rxns
TIGIT:CD112 Homogeneous Assay Kit	72030	384 rxns
TIGIT:CD155 Homogeneous Assay Kit	72029	384 rxns
PD-1:PD-L2 Homogeneous Assay Kit	72015	384 rxns

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com