



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## **Data Sheet**

### ***Blocking Buffer 2***

**Catalog #: 79728**

**Description:** This Blocking Buffer 2 is optimized for minimizing background in BPS Bioscience immunotherapy assay kits.

**Applications:** Suitable for use with enzymatic assays to study enzyme kinetics or screen for small molecular inhibitors in BPS Bioscience immunotherapy assay kits.

**Format:** Phosphate-Buffered solution

**Storage/Stability:** Store at 4°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 **1.858.202.1401** 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)