



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**

### ***5x SHP-2 Assay Buffer***

**Catalog #: 79626**

**Description:** The 5X SHP-2 Assay Buffer is optimized for use with the SHP-2 (Full-Length) Homogeneous Assay Kit (BPS Bioscience #79330) and the SHP-2 (catalytic domain) Homogeneous Assay Kit (BPS Bioscience #79317).

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

**Supplied As:** Tris-buffered solution

**Format:** Aqueous buffer solution

**Quantity:** 50 ml

**Storage/Stability:** Store at -20°C. Avoid freeze/thaw cycles. Stable 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)