



6042 Cornerstone Court West, Suite 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**

### ***Streptavidin-HRP*** Catalog # 79742

**DESCRIPTION:** Streptavidin-HRP is useful in BPS Bioscience assay kits.  
Note: This Streptavidin-HRP is *not* for use with BPS PARP assay kits.

**PROTOCOL:** Dilute Streptavidin-HRP in blocking buffer before use. See designated assay kit protocol for detailed dilution and assay instructions.

**APPLICATIONS:** Study enzyme kinetics, and screen small molecular inhibitors for drug discovery and HTS applications

**SUPPLIED AS:** Aqueous buffer solution

**STORAGE:** +4°C

**WARNING:** Avoid freeze/thaw cycles

**FORMAT:** Aqueous buffer solution

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)