



6042 Cornerstone Court West, Suite B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

Streptavidin-HRP (For PARP & Cytokine Assay Kits) Catalog # 80611

DESCRIPTION: Streptavidin-HRP is useful in select PARP and Cytokine BPS Bioscience assay kits.

PROTOCOL: Dilute Streptavidin-HRP 1:50 in blocking buffer before use. See designated assay kit protocol for detailed assay instructions.

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

SUPPLIED AS: Aqueous buffer solution

SIZE: 100 µl

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

Please visit our website at: www.bpsbioscience.com