



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

5X PTP Assay Buffer

Catalog #: 79716

DESCRIPTION: Extra assay buffer for use with the PTP Colorimetric assay kit (BPS Bioscience, Cat. #30019), or related PTP assay kits, for the study of protein tyrosine phosphatase activity.

APPLICATIONS: Ideal for detection of protein tyrosine phosphatase activity, and screening of small molecule inhibitors for drug discovery and HTS applications.

SUPPLIED AS: Aqueous buffer solution.

QUANTITY: 20 mL

STORAGE: Store in aliquots at -20°C.

STABILITY: At least 6 months at -20°C.

INSTRUCTIONS FOR USE: Please see appropriate assay kit protocol for detailed instructions.

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
10x PTP1B Colorimetric Substrate	79693	5 ml
PTP1B Colorimetric Assay Kit	30019	96 rxns

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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com