



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

10x PTP1B Colorimetric Substrate

Catalog #: 79693

Description: 10x PTP1B Colorimetric Substrate for use with BPS Biosciences' PTP1B Colorimetric Assay Kit (#30019).

Formulation: Aqueous solution of p-Nitrophenylphosphate

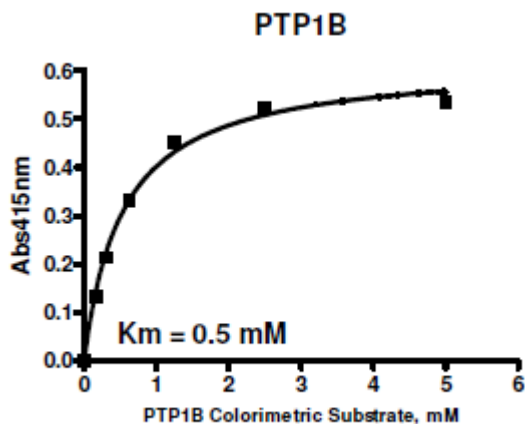
Concentration: 5 mM

Supplied as: Aqueous buffer solution

Size: 5 ml

Storage: -20°C

Example Assay Results:



References:

1. Protein-Tyrosine Phosphatase-1B Mediates Sleep Fragmentation-Induced Insulin Resistance and Visceral Adipose Tissue Inflammation in Mice (2017)

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