

Data Sheet

PTPmu

Human, recombinant, N-terminal GST tag
Catalog #: 30053
Lot#: 121108 **Conc.:** 0.35 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
 110 mM NaCl, 2.2 mM KCl, 16 mM
 glutathione, 20% glycerol, and 3 mM DTT.

Stability: >6 months at -80°C

References:

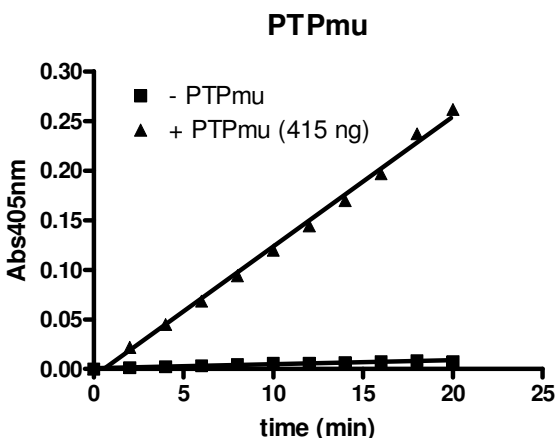
1. S. Brady-Kalnay, *et al.*, *J. Cell Biol* 1995, **130**: 977
2. T. Mourton, *et al.*, *J. Biol. Chem.* 2002, **276**: 14896

Description: PTPmu (protein tyrosine phosphatase mu, also known as RPTP μ), catalytic domain, (a.a. 896-1175), (GenBank Accession No. NM_001105244) with N-terminal GST tag, MW=56.2 kDa, expressed in an *E. coli* expression system.

Specific Activity: 691 pmole/min/ μ g.
Assay conditions: 20 mM HEPES, pH 7.2, 100 mM NaCl, 2 mM EDTA, 1 mM DTT, 10 mM pNPP and PTP-mu. Assay was done at 30°C for 15 min, and the reaction was monitored at 405 nm continuously. Specific activity was calculated based on $\epsilon_{\text{pNPP}^-} = 18,000 \text{ M}^{-1}\text{cm}^{-1}$

Application: Useful for the study of enzyme kinetics, regulation, to dephosphorylate target substrates and for screening inhibitors.

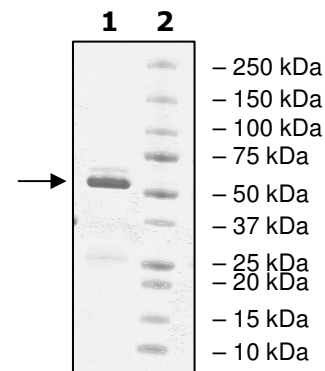
Quality Assurance



4-20% SDS-PAGE Coomassie staining

Lane 1:
3.36 μ g PTPmu
Lane 2:
Protein Marker

MW: 56.2 kDa
Purity: $\geq 81\%$



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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com