

Data Sheet

TC-PTP, GST-tag

Human, recombinant, N-terminal GST tag

Catalog #: 30013

Lot #: 150219

Conc.: 2.9 mg/ml

Formulated in: 25 mM Tris-HCl, pH 8.0, 75 mM NaCl, 0.05% Tween-20, 50% glycerol, 2 mM EDTA, 1 mM DTT, 10 mM glutathione.

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

1. N.F. Zander *et al.*, *Biochemistry* 1991, **30**: 6964.
2. D.E. Cool *et al.*, *Proc. Natl. Acad. Sci. USA* **86**: 5257.

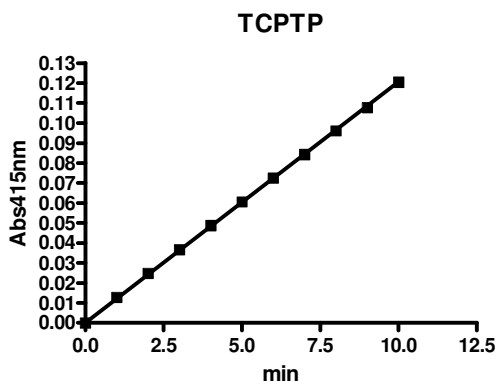
Description: Human recombinant TC-PTP, also known as PTPN2 (protein tyrosine phosphatase N2), catalytic domain, (a.a. 2-315), (Genbank accession number BC016727) with N-terminal GST tag, MW=62.3 kDa, expressed in an *E. coli* expression system.

Specific Activity: 1,000 pmole/min/ μg .
 Assay Conditions: Enzyme reaction (50 μl) containing 20 mM HEPES, pH 7.2, 100 mM NaCl, 2 mM EDTA, 1 mM DTT, 1 mM pNPP and PTPase was monitored at 415 nm continuously. Specific activity was calculated based on $\epsilon_{\text{pNP}^-} = 18,000 \text{ M}^{-1}\text{cm}^{-1}$

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

Quality Assurance

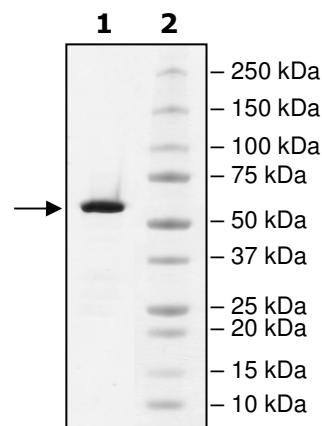
Specific Activity



4-20% SDS-PAGE
Coomassie staining

Lane 1:
 2 μg TC-PTP
Lane 2:
 Protein Marker

MW: 63.3 kDa
Purity: $\geq 90\%$



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