

Data Sheet

PEST(PTPN12)

Human recombinant, with N-terminal GST-tag

Catalog #: 30023

Lot #: 110215 Conc.: 1.2 mg/ml

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

Stability: >6 months at -80°C

References:

1. J.F. Cote, et al., *J Biol Chem.* 2002, **277**:
2973-86.
2. K. Horsch et al., *Mol Endocrinol.* 2001, **12**:
2182-96

Description:

Human recombinant PTP-PEST, also known as PTPN12, GenBank Accession # NM_002835), catalytic domain, (a.a. 2-300) with N-terminal GST tag, MW=61.4 kDa, expressed in *E. coli*.

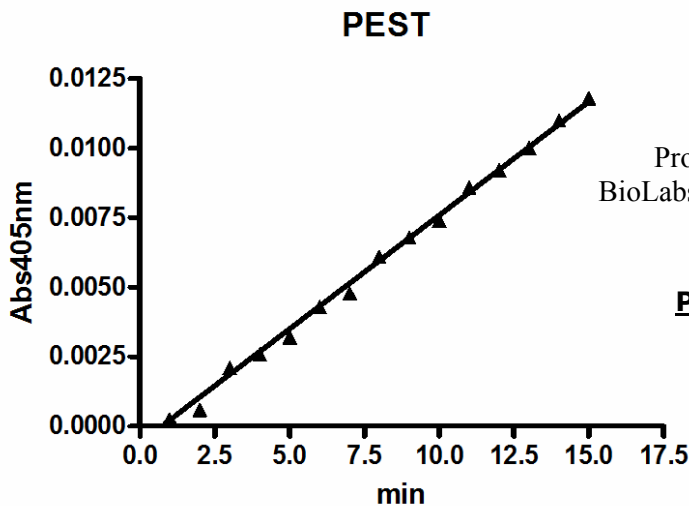
Specific Activity: 76 pmole/min/μg.

Assay condition: 20 mM HEPES, pH 7.2, 100 mM NaCl, 2 mM EDTA, 1 mM DTT, 1 mM pNPP and 2 ng/μl PEST. Assay was done at RT for 15 min, and the reaction was monitored at 405 nm continuously. Specific activity was calculated based on $\epsilon_{pNPP} = 18,000 \text{ M}^{-1}\text{cm}^{-1}$.

Application:

Useful for the study of enzyme kinetics, regulation, and to dephosphorylate target substrates.

Quality Assurance



10% SDS-PAGE Coomassie staining

Lane 1:

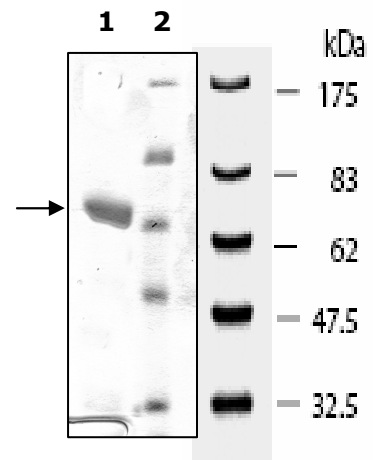
4 μg PEST

Lane 2:

Protein Marker
BioLabs (#P7708L)

MW: 61 kDa

Purity: >80%



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