

Data Sheet

PTP IA2

Human, recombinant, N-terminal GST-tag

Catalog #: 30054

Lot #: 1007

Conc.: 0.4 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, and 10 mM glutathione.

Stability: >6 months at -80°C

References:

1. Ort, T., *et al.*, *EMBO J.* 2001, **20**: 4013-23
2. Li, Z., *et al.*, *J Histochem Cytochem.* 2000, **48**:761-7

Description:

Protein tyrosine phosphatase. (PTP-IA2, ICA-512), Catalytic domain, (a.a. 693-979), (Gene bank accession no. 002846), with N-terminal GST tag, MW=58.8 kDa, expressed in an *E. coli* expression system.

Application:

Suitable for use as an immunogen, in binding assays or as an antigen in autoimmune assays.

Quality Assurance

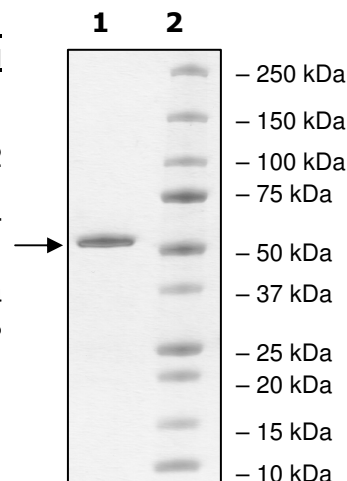
4-20% SDS-PAGE Coomassie staining

Lane 1:
2 µg PTP-IA2

Lane 2:
Protein Marker

MW: 58.8 kDa

Purity: ≥95%



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 **1.858.259.5734**

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

Please visit our website at: www.bpsbioscience.com