

Data Sheet

SHP-2, His-Tag (Full Length)

Human, Recombinant, N-terminal His-tag
Catalog #: 79018
Lot #: 170328-G2 **Conc.:** 0.41 mg/ml

Formulated in: 40mM Tris pH8.0, 110mM NaCl, 2.2mM KCl, 0.04% Tween-20 and 20% glycerol.

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Protein may be diluted to $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at -80°C .

References:

1. B.G. Neel *et al.* *Trends Biochem. Sci.* (2003) **28**: 284
2. J.V. Ravetch & L.L. Lanier *Science* (2000) **290**: 84

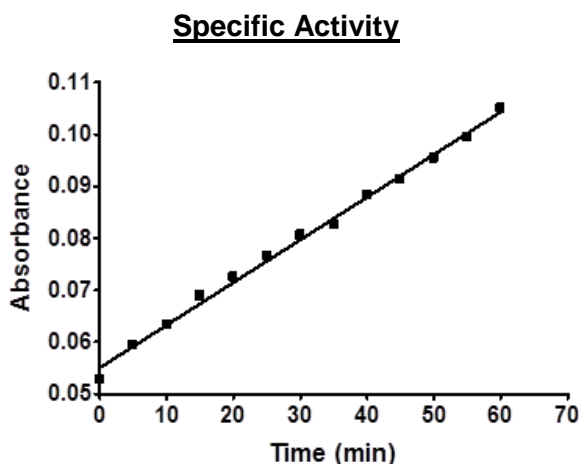
Description: Human recombinant SHP-2, also known as PTPN11 (protein tyrosine phosphatase N11), a.a. 2-593 (end), full length, with N-terminal His-tag, Genbank accession number NM_002834, MW=70 kDa, expressed in an *E. coli* expression system.

Specific Activity: $\geq 50 \text{ pmol/min}/\mu\text{g}$.

Assay conditions: 50mM HEPES, pH7.4, 100mM NaCl, 2mM EDTA, 3mM DTT, 5mM PNPP substrate, 30 min at 30°C . Absorbance is measured at 405nm.

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

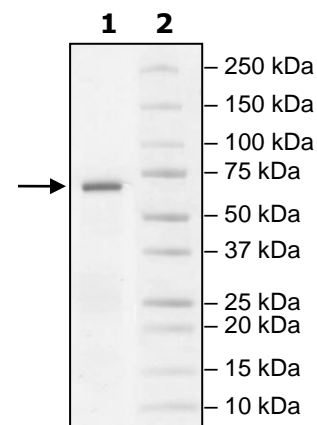


**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
2 μg SHP2

Lane 2:
Protein Marker

MW: 70 kDa
Purity: $\geq 85\%$



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