

Data Sheet

Jak2 (JH1, JH2 Domain), His-GST-Tag

Human Recombinant, N-terminal His-GST-tag
Catalog #: 40451
Lot #: 190115-2 **Conc.:** 0.15 mg/ml

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

Stability: At least 6 months at -80°C .
 Avoid freeze/thaw cycles.

References:

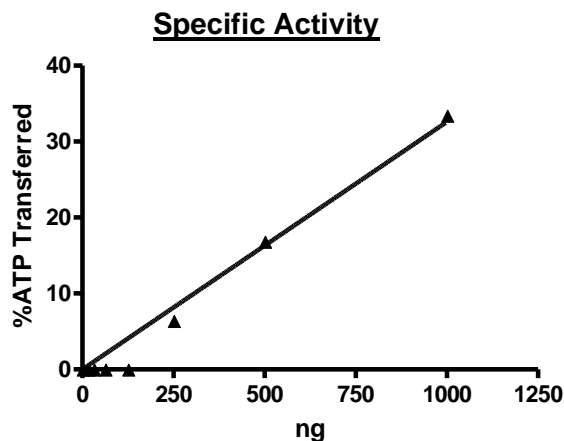
1. Tono, C., *et al.*, *Leukemia* **19(10)**: 1843-1844 (2005).
2. Goertler, P.S., *et al.*, *Blood* **106(8)**: 2862-2864 (2005).

Description: Human Jak2, also known as Janus Kinase 2, Tyrosine-Protein Kinase JAK2, and THCYT3, GenBank Accession No. NM_004972, containing the catalytic and the pseudo kinase domains, a.a. 532-1132 (end), with N-terminal His-GST-tag, expressed in a Baculovirus infected Sf9 cell expression system. MW=98 kDa.

Specific Activity: ≥ 12 pmol/min/ μg
 Assay Conditions: 40 mM Tris-HCL pH 7.4, 20 mM MgCl_2 , 0.1 mg/mL BSA and 0.2 mM DTT using 0.2 mg/ml Poly-(Glu4:Tyr) substrate and 20 μM of ATP. Reaction was done at 30°C for 45 min. Amount of ATP transferred was calculated using Kinase-Glo reagent from Promega.

Application: Useful for studying enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

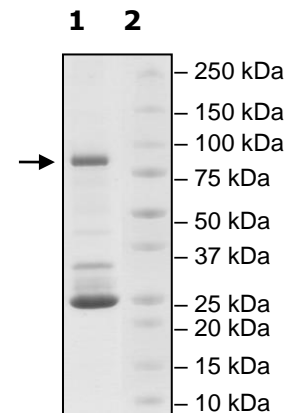


**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
JAK2
Lane 2:
 Protein Marker

Purity: $\geq 25\%$

MW: 98 kDa



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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com