

Data Sheet

Jak2 (V617F), GST-Tag, Avi-Tag, Biotin-Labeled (Dephosphorylated)

Human, Recombinant, N-terminal GST-Tag
 C-terminal Avi-Tag, Biotin Labeled
Catalog #: 79115
Lot #: 200701 **Conc.:** 0.06 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-
 20, 16 mM Glutathione, 3mM DTT, 20%
 glycerol.

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

References:

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

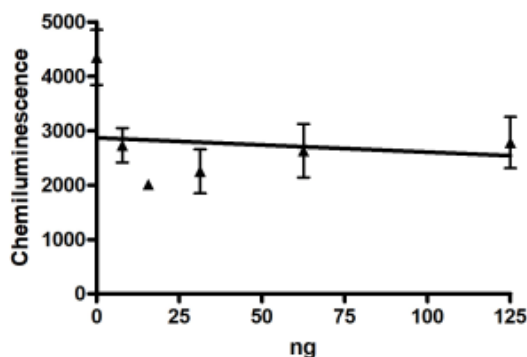
Description: Human Jak2 containing the catalytic and the pseudo kinase domain, also known as Janus kinase-2, Just Another Kinase 2, Jak-2-JH1 and JH2, GenBank Accession No. NM_004972, a.a. 513-1132(end), with V617F mutation; with N-terminal GST-Tag and C-terminal Avi-Tag, co-expressed with PTP1B for dephosphorylation in HEK293 expression system. This protein is enzymatically Biotin-labeled using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%. MW=100 kDa.

Assay Conditions: Assay was performed according to the TYR Phosphorylation Assay Kit (BPS Bioscience, #79502), with both phosphorylated JAK2 (standard) and JAK2 titrated from 500 ng to 0 ng.

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

Quality Assurance

JAK2 Phosphorylation Detection

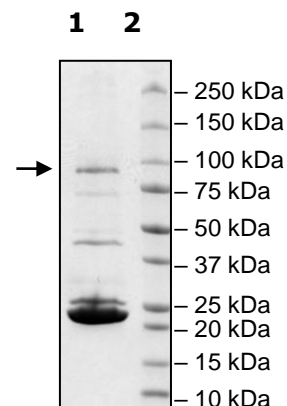


4-20% SDS-PAGE Coomassie staining

Lane 1:
JAK2
Lane 2:
 Protein Marker

Purity: ≥8%

MW: 100 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 **1.858.259.5734**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com