

Data Sheet

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

Human, Recombinant, N-terminal GST-Tag
C-terminal Avi-Tag, Biotin Labeled

Catalog #: 79115

Lot #: 171213

Conc.: 0.2 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 0.04% Tween-
20, 20mM GSH and 20% glycerol.

Stability: At least 6 months at -80°C . *Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).*

References:

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

Description:

Human Jak2 (JH1 catalytic domain and JH2 pseudokinase domain) with V617F mutation, GenBank Accession No. NM_004972, a.a. 513-1132 (end), with N-terminal GST tag, C-terminal Avi-Tag, expressed in a HEK293 cell expression system and enzymatically Biotin-labeled using AviTag™ technology. MW=100 kDa. No phosphorylation was detected by ELISA.

Quality Assurance

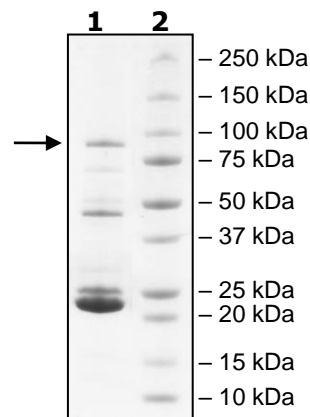
4-20% SDS-PAGE Coomassie staining

Lane 1:
0.5µg JAK2(V617F)

Lane 2:
Protein Marker

MW: 99.5 kDa

Purity: ≥10%



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** or **1.858.481.8694**, Fax **1.858.259.5734**

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