

Data Sheet

Jak2, FLAG-Avi-Tag, Biotin-Labeled

Human, Recombinant, C-terminal
FLAG-Avi-Tag, Biotin Labeled

Catalog #: 79493

Lot #: 190913

Conc.: 0.29 mg/ml

***Concentration listed is for purity %
adjusted and biotin-labeled target protein**

Description: Full length human Jak2, also known as Janus kinase-2, Just Another Kinase 2, JAK2-JH2, GenBank Accession No. NM_004972, a.a. 1-1132(end), with C-terminal FLAG-Avi-Tag, co-expressed with PTP1B in HEK293 expression system. This protein is enzymatically Biotin-labeled using Avi-Tag™ technology. MW=134 kDa. Phosphorylation is confirmed to be <15%. Biotinylation is confirmed to be ≥90%.

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 0.04% Tween-
20, 100 µg/ml FLAG peptide and 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. Goerttler, P.S., *et al.*, *Blood*. 2005;
106(8): 2862-2864.

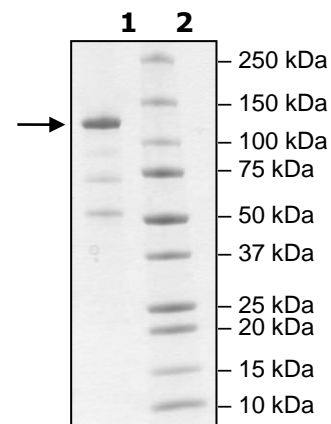
Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
2µg JAK2
Lane 2:
Protein Marker

MW: 134 kDa

Purity: ≥75%



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