

Data Sheet

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

Human, Recombinant,
C-terminal His-Avi-Tag, Biotin Labeled
Catalog #: 79075
Lot #: 180514 **Conc.:** 0.13 mg/ml

Description:

Human Jak2, also known as Janus kinase-2, Just Another Kinase 2, with V617F mutation, GenBank Accession No. NM_004972, a.a. 513-827, with C-terminal His-Avi-Tag, expressed in an Sf9 expression system. This protein is enzymatically Biotin-labeled using Avi-Tag™ technology. MW=39 kDa.

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 500 mM Imidazole 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%).

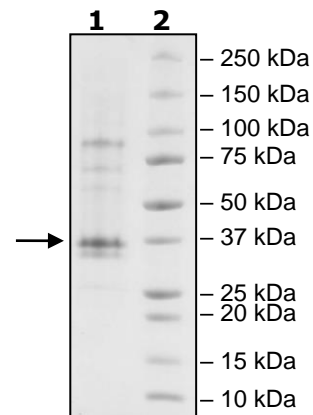
References:

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

Quality Assurance

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

Lane 1:
2 μg JAK2(V617F)
(513-827-
His-Avi-Biotin)
Lane 2:
Protein Marker
MW: 39 kDa
Purity: $\geq 58\%$



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