

Data Sheet

Jak2 (JH2 Domain) (E596A, V617F, W659A, W777A, F794H), His-Tag

Human, Recombinant, C-terminal His-tag

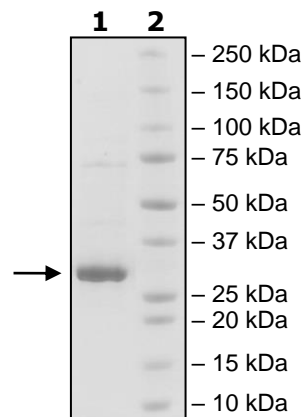
Catalog #: 100122**Lot #:** 180725-G1 **Conc.:** 0.19 mg/ml**Formulated in:** 20 mM Tris pH 8.5, 0.5 mM TCEP and 10% glycerol**Stability:** At least 6 months at -80°C . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).**References:**

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 (JH2 pseudokinase domain), also known as Janus kinase-2, and Just Another Kinase 2, with E596A, V617F, W659A, W777A, and F794H mutations. GenBank Accession No. NM_004972, a.a. 535-812, with C-terminal His-Tag, MW=33 kDa, expressed in a Sf9 expression system.

Quality Assurance

**4-20% SDS-PAGE
Coomassie staining****Lane 1:**
2 μg JAK2 (E596A,
V617F,W659A,
W777A,F794H)**Lane 2:**
Protein Marker
MW: 33 kDa
Purity: $\geq 90\%$ 

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