

Data Sheet

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

Human, Recombinant, C-terminal His-tag

Catalog #: 100123

Lot #: 180829

Conc.: 2.0 mg/ml

Formulated in: 20 mM Tris-HCl, pH 8.5,
10% glycerol, 0.5mM TCEP

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. Goertler, 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 W659A, W777A, and F794H mutations. GenBank Accession No. NM_004972, a.a. 536-812, with C-terminal His-Tag, MW=33 kDa, expressed in a Sf9 expression system.

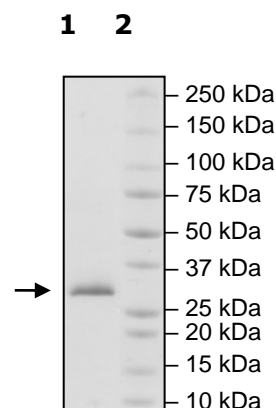
Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
JAK2-3 mutations

Lane 2:
Protein Marker
Purity: >90%

MW: 33 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** 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