

Data Sheet

Jak2 (W659A, W777A, F794H), His-Avi-Tag, Biotin

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

Catalog #: 79463

Lot #: 180323

Conc.: 1.86 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 0.04% Tween-
20, 140 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.

Description:

Human Jak2 (JH1 catalytic domain and
JH2 pseudokinase domain) with W659A,
W777A, and F794H mutations, GenBank
Accession No. NM_004972, a.a. 536-
812, with C-terminal His-Avi-Tag,
expressed in an Sf9 cell expression
system and enzymatically Biotin-labeled
using AviTag™ technology. MW=35
kDa.

Quality Assurance

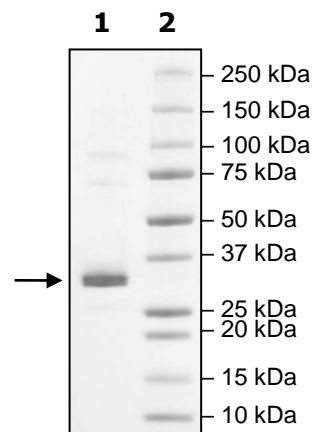
4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg JAK2(W659A,
W777A, F794H)

Lane 2:
Protein Marker

MW: 35 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