



6042 Cornerstone Court West, Suite B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

Jak2, His-tag

Human, Recombinant, N-terminal His-tag

Catalog #: 40450

Lot #: 180411-G **Conc.:** 0.25 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 3 mM DTT,
0.04% Tween-20 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 (GenBank Accession No.
NM_004972), (a.a. 808-end) with N-
terminal His tag, MW=43 kDa,
expressed in Baculovirus expression
system.

Specific Activity: 124 pmol/min/ μg

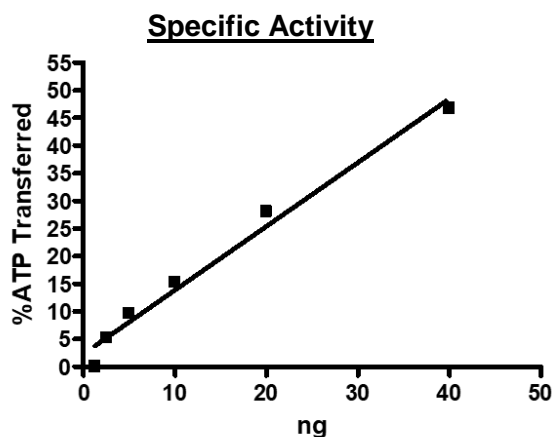
Assay Conditions: Assay was done in
Kinase buffer containing 1 mM DTT
using Poly-(Glu4:Tyr) substrate (0.2
mg/ml) and 20 μM ATP. Reaction was
done at 30°C for 45 min.

Amount of ATP transferred was
calculated using Kinase-Glo reagent
(Promega)

Application:

Useful for the study of enzyme kinetics,
screening inhibitors, and selectivity
profiling.

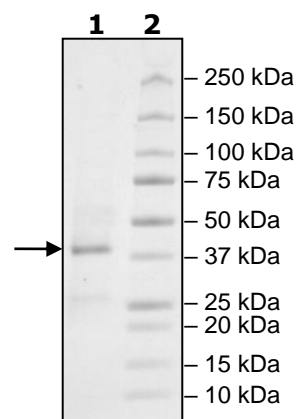
Quality Assurance



4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg JAK2
Lane 2:
Protein Marker

MW: 43 kDa
Purity: $\geq 71\%$



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