

Data Sheet

Jak3, His-tag

Human, recombinant, N-terminal 10x His-Tag
Catalog #: 40452
Lot #: 150921-A2 **Conc.:** 0.79 mg/ml

Formulated in: 40 mM Tris pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 200 mM imidazole, 3 mM DTT and 20% glycerol

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles.

References:

1. Kim S. *et al.*, *Cell. Signal.* **21 (6)**, 892-898 (2009).
2. Mullighan C.G. *et al.*, *Proc. Natl. Acad. Sci. U.S.A.* **106 (23)**, 9414-9418 (2009).

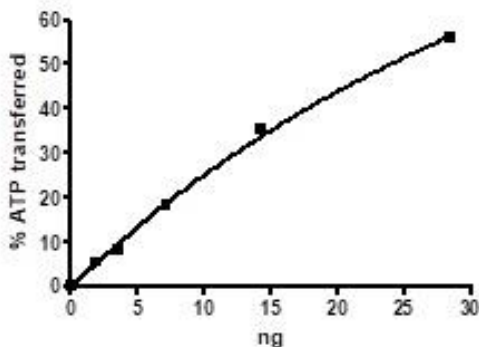
Description: Human Jak3 (Janus kinase 3), also known as LJAK, GenBank Accession No. NM_000215, a.a. 811-1124(end) with N-terminal 10x His-Tag, expressed in Baculovirus-infected Sf9 cell expression system. MW = 37 kDa.

Specific Activity: 500 pmol/min/ μg .
Assay conditions: 40 mM Tris-HCL pH 7.4, 20 mM MgCl₂, 0.1 mg/mL BSA and 0.2 mM DTT using 0.2 mg/ml Poly-(Glu4:Tyr) substrate and 20 mM ATP. Reaction was done at 30°C for 45 min. Amount of ATP transferred was calculated using Kinase-Glo reagent from Promega.

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

Quality Assurance

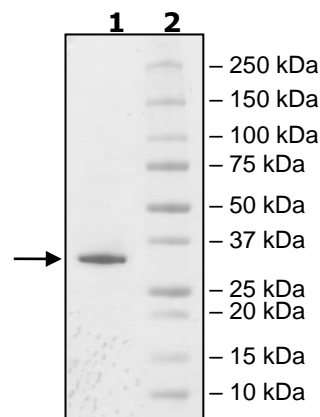
Specific Activity



4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg JAK3
Lane 2:
Protein Marker

MW: 37 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** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com