

## Data Sheet

### Jak2 (JH1, JH2 domain), GST-Tag, Avi-Tag, Biotin

Human, Recombinant, N-terminal GST-Tag,  
C-terminal Avi-Tag, Biotin Labeled

**Catalog #:** 79495

**Lot #:** 180507

**Conc.:** 0.023 mg/ml

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

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . *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.

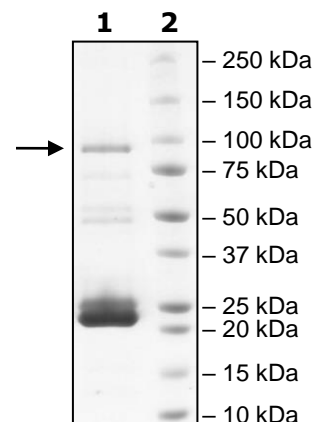
#### Description:

Human Jak2 containing the catalytic and the pseudo kinase domain (a.a. 513-1132), also known as Janus kinase-2, Just Another Kinase 2, Jak-2-JH1, JH2, GenBank Accession No. NM\_004972, with N-terminal GST-Tag and C-terminal Avi-Tag™ co-expressed with PTP1B in HEK293 expression system. This protein is enzymatically Biotin-labeled using AviTag™ technology. Percentage of phosphorylation  $\leq 10\%$ . MW=100 kDa.

## Quality Assurance

### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
1 $\mu\text{g}$  JAK2  
**Lane 2:**  
Protein Marker  
**MW:** 100 kDa  
**Purity:**  $\geq 10\%$



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](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)