

## Data Sheet

### Jak1, GST-Tag

Human, recombinant, N-terminal GST-Tag

**Catalog #:** 40449

**Lot#:** 190919-3

**Conc.:** 0.20 mg/ml

**Formulated in:** 25 mM Tris-HCl pH 8.0, 427 mM NaCl, 1.3 mM KCl, , 0.02% Tween 20, 3 mM DTT, 14 mM glutathione and 50% glycerol.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles.

**References:**

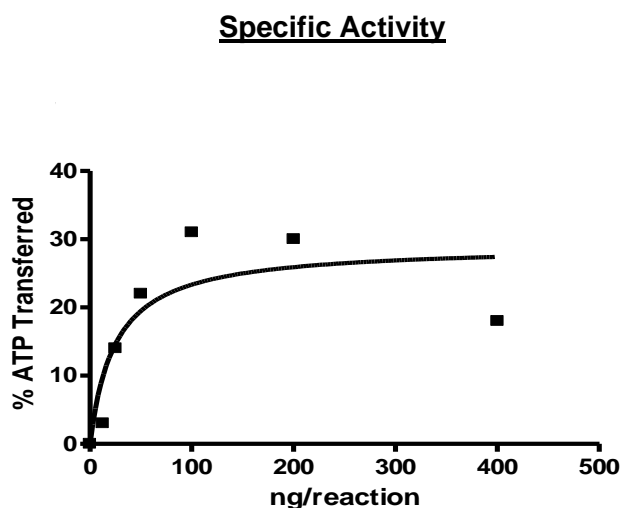
1. Rodig, S.J., *et al.* (1998). *Cell* **93 (3)**: 373–8.
2. Haan, C., *et al.* (2001). *J. Biol. Chem.* **276 (40)**: 37451–8.
3. Miyazaki, T., *et al.* (1994). *Science* **266 (5187)**: 1045–7.

**Description:** Human Jak1 also known as Janus kinase 1 and JTK3, GenBank Accession No. NM\_002227, a.a. 866-1154(end) with an N-terminal GST-Tag, expressed in Sf9 cells via a baculovirus expression system, MW= 60 kDa.

**Specific Activity:**  $\geq 80$  pmole/min/ $\mu\text{g}$ . Assay was done in Kinase buffer containing 2 mM DTT using IRS-1 derived peptide substrate (0.1 mg/ml) and 20  $\mu\text{M}$  ATP. Reaction was done at  $30^{\circ}\text{C}$  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

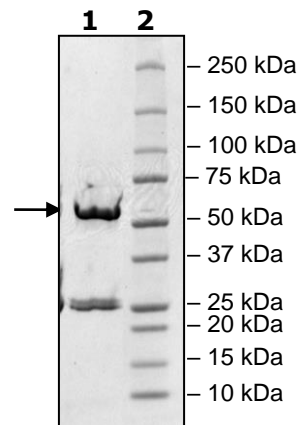


**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
4  $\mu\text{g}$  JAK1  
**Lane 2:**  
Protein Marker

**MW:** 60 kDa

**Purity:**  $\geq 63\%$



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](mailto:info@bpsbioscience.com)

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