

## Data Sheet

### Tyk2, His-Tag Recombinant

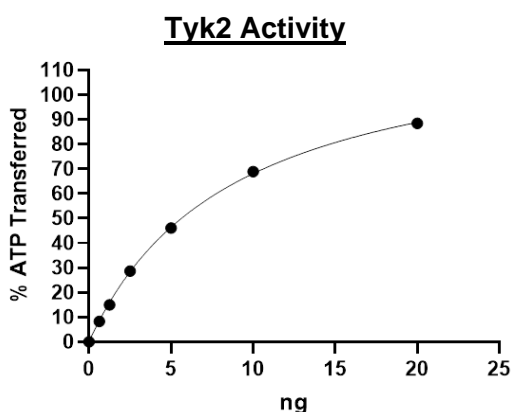
Human, recombinant, N-terminal His-tag

**Catalog #:** 40285**Lot#:** 110830      **Conc.:** 0.18 mg/ml**Formulated in:** 45 mM Tris, pH 8, 125 mM NaCl, 2.4 mM KCl 0.045% Tween-20, 100 mM imidazole and 10% glycerol.**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles.**References:**

1. Firmbach-Kraft I, *et al.* (1990). *Oncogene* **5 (9)**: 1329–36.
2. Partanen J, *et al.* (1991) *Proc. Natl. Acad. Sci. U.S.A.* **87 (22)**: 8913–7.
3. Domanski P, *et al.* (1995). *FEBS Lett.* **374 (3)**: 317–22.

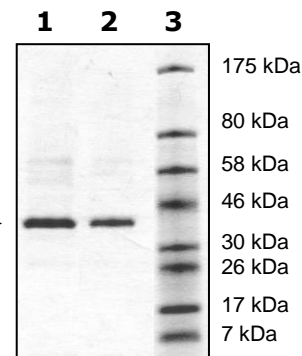
**Description:** Human Tyrosine Kinase 2 (Tyk2, also known as JTK1) GenBank Accession No. NM\_003331, a.a. 871-1187(end), with N-terminal His-tag, expressed in a Baculovirus-infected Sf9 cell expression system. MW=38 kDa.**Assay Conditions:** Assay was done according to TYK2 (Tyrosine Kinase 2) Assay Kit (BPS Bioscience #79527) with various amounts of Tyk2.**Application:** Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Assurance

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

**Lane 1 & 2:**  
4 & 2  $\mu\text{g}$  Tyk2  
**Lane 3:**  
Protein Marker  
BioLabs (#P7708L)

**MW:** 38 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: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)