

Data Sheet

Tyk2(JTK1)

Human, recombinant, N-terminal His-tag

Catalog #: 40285

Lot#: 100930 Conc.: 0.6 mg/ml

Formulated in:

50 mM Tris, pH 8, 130 mM NaCl, 2.4 mM KCl
0.04% Tween-20, 250 mM imidazole and 15%
glycerol.

Stability:

>6 months at -80°C

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.

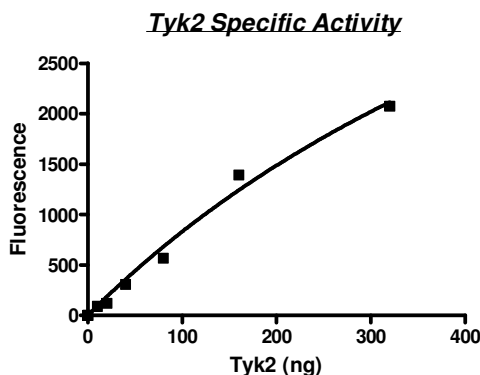
Description:

Human Tyk2 (Tyrosine Kinase 2, also known as JTK1), GenBank Accession No. NM_003331, a.a. 871-1187(end), with N-terminal His-tag, MW=38 kDa, expressed in a Baculovirus-infected Sf9 cell expression system.

Assay conditions: Add 50 µl reaction mix (Tyk2 diluted in 1x kinase buffer supplemented with 0.2 mM ATP) to wells coated with poly(Glu-Tyr, 4:1) substrate. Incubate for 1h at RT followed by addition of Europium-labeled anti-phospho-Tyr antibody and incubate for another 1h at room temperature. Time resolved fluorescence measured using development solution at $\lambda_{exc}=330\pm 40$ nm and $\lambda_{em}=620\pm 10$ nm.

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

Quality Assurance



10% SDS-PAGE Coomassie staining

Lane 1:

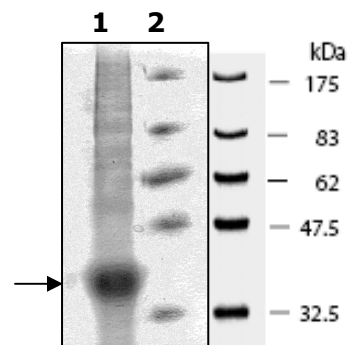
8 µg Tyk2

Lane 2:

Protein Marker
BioLabs (#P7708L)

MW: 38 kDa

Purity: >55%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com