

Data Sheet

TAK1-TAB1

Human recombinant, with N-terminal GST-tag

Catalog #: 40279

Lot #: 120822-G2 Conc.: 0.58 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 3 mM DTT and 20% glycerol.

Stability: 12 months at -70°C

References:

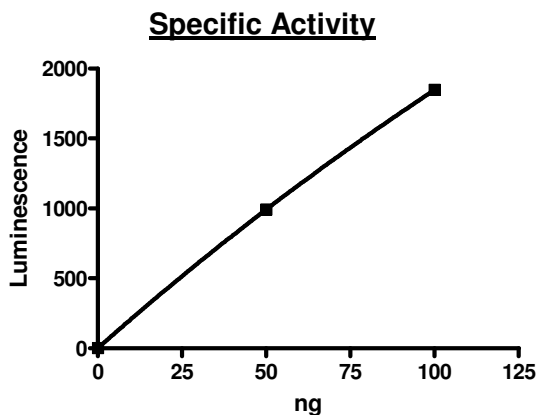
1. Sakurai, H., *et al.* (2000). *FEBS Lett. (Netherlands)* **474 (2-3)**: 141-5.
2. Liu, H.-H., *et al.* (2006). *Proc. Natl. Acad. Sci. USA* **103**: 11677-11682.
3. Yamaguchi, K., *et al.* (1995). *Science* **270**: 2008-2011.

Description: Recombinant human TAK1, also known as MAP3K7 (Mitogen-activated protein kinase kinase kinase 7), Genbank Accession #NM_003188, a.a. 1-303, and human TAB1, Genbank Accession #NM_006116, a.a. 437-end, linked by a small peptide (DFGGGGG), and with an N-terminal GST-tag, MW = 69 kDa, expressed in Sf9 cells via a baculovirus expression system.

Assay Conditions: Assay was performed in a kinase buffer containing 0.2 mM DTT using MBP substrate (0.1 mg/ml) and 100 mM ATP. Reaction was done at 30°C for 35 min. Amount of ATP transferred was calculated using ADP-Glo® reagent (Promega).

Application: Useful for studying enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

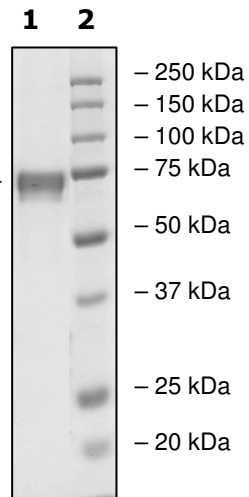


4-20% SDS-PAGE Coomassie staining

Lane 1:
2.5 μg TAK1-TAB1

Lane 2:
Protein Markers

MW: 69 kDa
Purity: $\geq 93\%$



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