

Data Sheet

JNK1

Mouse, recombinant, N-terminal GST tag

Catalog #: 40071

Lot #: 111111 **Conc.:** 0.1 mg/ml

Formulated in: 50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 0.1 mM EDTA, 0.1 mM EGTA, 0.25 mM DTT, 0.1 mM PMSF, and 25% glycerol.

Stability: >12 months at -80°C

References:

1. Derijard, B. *et al.*, Cell. 1994 Mar 25; **76(6)**: 1025-1037.
2. Kolodziejczyk, S M. *et al.*, Curr. Biol. 2001 Aug 21; **11(16)**: 1278-1282.

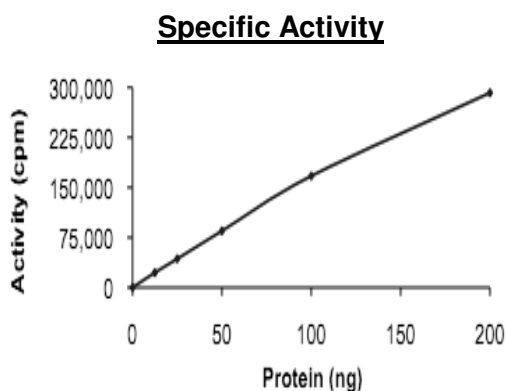
Description: Mouse JNK1, also known as SAPK1 or PRKM8 (GenBank Accession No. NM_016700), full length, with N-terminal GST tag, MW= \sim 70 kDa, expressed in Sf9 insect cells via a baculovirus expression system.

Specific Activity: 93 pmol/min/ μ g

Assay Buffer: 5 mM MOPS, 2.5 mM β -glycerophosphate, 5 mM MgCl₂, 1 mM EGTA, 0.4 mM EDTA, 0.05 mM DTT, 2 mM ATP. Incubate JNK1 with ATF2 substrate (0.1 mg/ml) and [³³P]-ATP at 30°C for 15 minutes, then spot reaction on phosphocellulose paper, fix in 1% phosphoric acid, and assay with a scintillation counter.

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

Quality Assurance

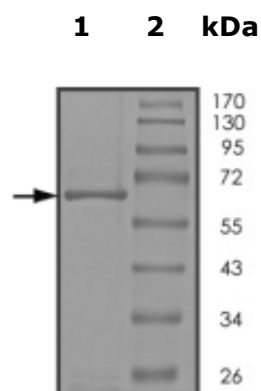


10% SDS-PAGE
Coomassie staining

Lane 1:
JNK1

Lane 2:
Protein Marker

MW: \sim 70 kDa
Purity: \geq 95%



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