

Data Sheet

JNK1- β 1(K55M)

Human, recombinant, N-terminal GST tag

Catalog #: 40871

Lot #: 100119 **Conc.:** 2.4 mg/ml

Formulated in: 45 mM Tris-HCl, pH 8.0,
124 mM NaCl, 2.4mM KCl, 10% glycerol,
and 3 mM DTT.

Stability: >12 months at -80°C

References:

1. Derijard, B. et al., *Cell*. **76(6)**:1025-37 (1994).
2. Kolodziejczyk, S M. et al., *Curr. Biol.* **11(16)**:1278-82 (2001).

Description:

Human JNK1- β 1/MAPK8 (K55M) (GenBank Accession No. NM_139046), (a.a. 2-383) with N-terminal GST tag, MW= \sim 70 kDa, expressed in *E. coli*.

Application:

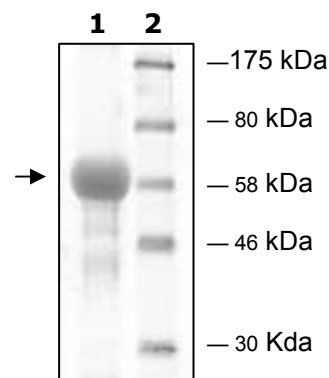
Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

**10% SDS-PAGE
Coomassie staining**

Lane 1:
14 μ g JNK1- β 1

Lane 2:
Protein Marker
BioLabs (#P7708L)



MW: \sim 70 kDa

Purity: \geq 85%

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