

ProQinase™ JNK1 K55R/K56R

mitogen-activated protein kinase 8

Recombinant Protein Kinase Substrate

HGNC Symbol: MAPK8

Synonyms: PRKM8, SAPK1

Product No.: 0524-0000-1

Lot: 006

Description: Human JNK1, full length, amino acids M₁-Q₃₈₄ (as in [NCBI/Protein](#) entry NP_002741.1), inactivated by K55R and K56R mutations, N-terminal HIS₆ fusion protein with a TEV cleavage site, expressed in E.coli

Theoretical MW_{Fusion Protein}: 46,466 Da

Expression host: E.coli

Purification: Immobilized Metal Affinity Chromatography

Storage buffer: 50 mM HEPES pH 7.5, 100 mM NaCl, 1 mM DTT, 10% glycerol

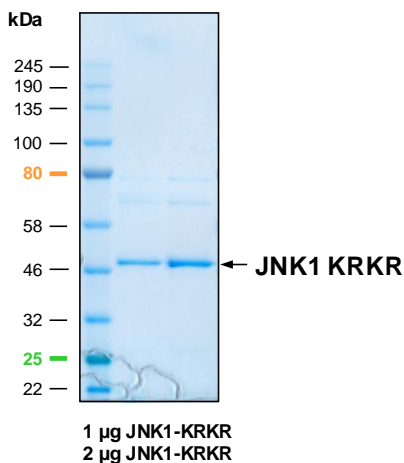
Storage temperature: -80°C

For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

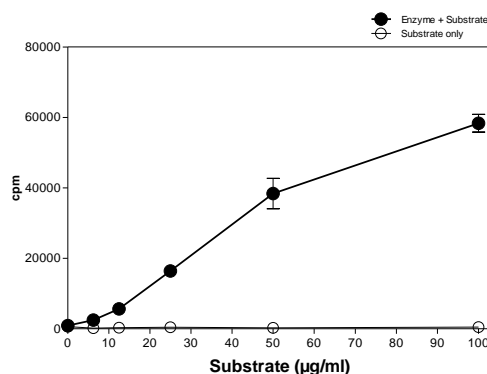
Protein concentration: 1.758 µg/µl

(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Coomassie stain



Phosphorylation of JNK1 K55R/K56R by the kinase MKK4 (Radiometric filter binding assay):



Assay conditions:

70 mM HEPES-NaOH, pH 7.5
3 mM MgCl₂
3 mM MnCl₂
3 µM Na-orthovanadate
1.2 mM DTT
50 µg/ml PEG_{20,000}
ATP: 1 µM
Substrate: variable concentration
Kinase: 1 µg/ml
MSFC membrane (Millipore)

Field of application:

JNK1 K55R/K56R has been validated for use in radiometric in-vitro kinase activity assays. It was not validated for use in ATP-consumption based kinase activity assays.

ProQinase™ JNK1 K55R/K56R

Product No.: 0524-0000-1

JNK1 K55R/K56R Recombinant Fusion Protein Amino Acid Sequence								
1	MHHHHH	AMT	ENLYFQG	AMS	RSKRDN	FYS	VEIGDSTFTV LKRYQNLKPI GSGAQGIVCA	60
61	AYDAILERNV	AI	RRLSRPFQ	NQTHAKRAYR	ELVLMKCVNH	KNIIGLLNVF	TPQKSLEEFQ	120
121	DVYIVMELMD	ANLCQVIQME	LDHERMSYLL	YQMLCGIKHL	HSAGIIHRDL	KPSNIVVKS	D	180
181	CTLKILDFGL	ARTAGTSFMM	TPYVVTRYR	APEVILGMGY	KENVDLWSVG	CIMGEMVCHK		240
241	ILFPGRDYID	QWNKVIEQLG	TPCPEFMKKL	OPTVVRTYVEN	RPKYAGYSFE	KLFPDVLFP	A	300
301	DSEHNKLGAS	QARDLLSKML	VIDASKRISV	DEALQHPYIN	VWYDPSEAEA	PPPKIPDKQL		360
361	DEREHTIEEW	KELIYKEVMD	LEERTKNGVI	RGQPSPLAQV	QQ			420

Red: 6HIS-tag Green: TEV site blue: JNK1 K55R/K56R boxed: K55R/K56R point mutations

JNK1 wt1 Amino Acid Sequence								
1	MSRSKRDN	FYS	VEIGDSTF	TVLKRYQNLK	PIGSGAQGIV	CAAYDAILER	NVAIKKLSRP	60
61	FQNQTHAKRA	YRELVLMKCV	NHKNIIGLLN	VFTQKSLEE	FQDVYIVMEL	MDANLCQVIQ		120
121	MELDHERMSY	LLYQMLCGIK	HLHSAGIIHR	DLKPSNIVVK	SDCTLKILDF	GLARTAGTSF		180
181	MMPYVVTRY	YRAPEVILGM	GYKENVDLWS	VGCIMGEMVC	HKILFPGRDY	IDQWNKVIEQ		240
241	LGTPCPEFMK	KLQPTVRYV	ENRPKYAGYS	FEKLPDVLV	PADSEHNKLG	ASQARDLLSK		300
301	MLVIDASKRI	SVDEALQHPY	INVWYDPSEA	EAPPPKIPDK	QLDEREHTIE	EWKELIYKEV		360
361	MDLEERTKNG	VIRGQPSPLA	QVQQ					420

blue: JNK1 sequence expressed in fusionprotein Red: variant in fusionprotein

¹NCBI/Protein accession number NP_002741.1