

ProQinase™ JNK1 (non-activated)

mitogen-activated protein kinase 8

Recombinant Human Protein Kinase

HGNC Symbol: MAPK8

Synonyms: PRKM8, SAPK1, JNK

Product No.: 0458-0000-2

Lot: 007

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

Theoretical MW_{Fusion Protein}: 46,410 Da

Expression host: E.coli

Purification: Immobilized Metal Affinity Chromatography

Activation: This kinase was not activated by special procedures

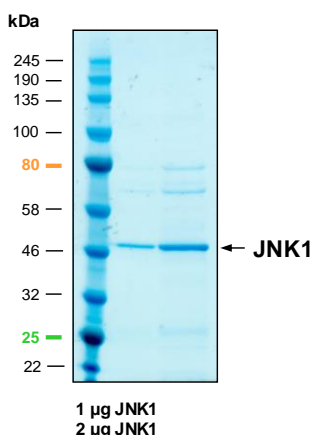
Storage buffer: 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 20 % glycerol

Storage temperature: -80°C

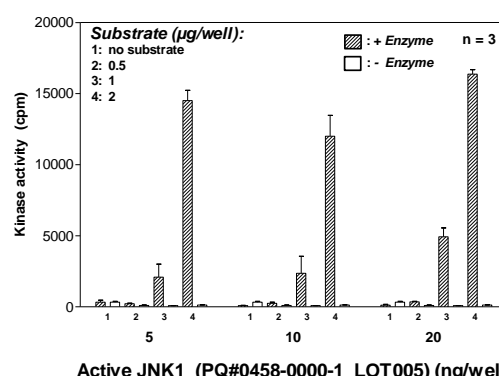
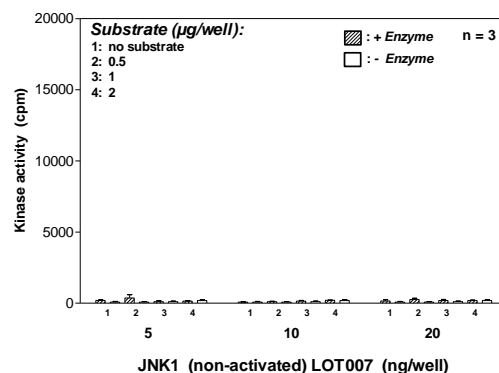
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 0.732 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

JNK1 Lot 007:
Coomassie stain



Kinase activity JNK1 (non-activated) vs active JNK1:



- Assay conditions:
60 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}
1 µM ATP
Substrate: ATF2 (variable concentration)
Kinase: JNK1 activated/non-activated
- Assay: ³³PanQinase® Assay

ProQinase™ JNK1 (non-activated)

Product No.: 0458-0000-2

HIS-JNK1 Recombinant Fusion Protein Amino Acid Sequence													
1	MHHHHH	AMT	ENLYFQG	AMS	RSKRDN	NFY	VEIGD	STFTV	LKRYQ	NLKPI	GSGAQ	GIVCA	60
61	AYDAI	LERNV	AIKKL	SRPFQ	NQTHA	KRAYR	ELVLM	KCVNH	KNIIG	LLNVF	TPQKS	LLEEFQ	120
121	DVYIV	MELMD	ANLCQ	VIQME	LDHER	MSYLL	YQMLC	GIKHL	HSAGI	IHRDL	KPSNI	VVKS	180
181	CTLKI	LDFGL	ARTAG	TSFMM	TPYVV	TRYR	APEVI	LGMGY	KENVD	LWSVG	CIMGE	MVCHK	240
241	ILFPG	RDRYD	QWNKV	IEQLG	TPCPE	FMKKL	QPTVR	TYVEN	RPKYA	GYSFE	KLFPD	VLFPA	300
301	DSEHN	KLKAS	QARDL	LSKML	VIDASK	RISV	DEALQ	HPYIN	VWYDP	SEAEA	PPPKI	PKQL	360
361	DEREH	TIEEW	KELIY	KEVMD	LEERT	KNGVI	RGQPS	PLAQV	QQ				420

Red: HIS6-tag Green: TEV cleavage site blue: JNK1

JNK1 wt ¹ Amino Acid Sequence													
1	MSRSKR	DNF	YSVEI	GDSTF	TVLKRY	QNLK	PIGSGA	QGIV	CAAYDA	ILER	NVAIK	LSRP	60
61	FQNQTH	AKRA	YRELV	LKCV	NHKNII	GLLN	VFTPK	SLEE	FQDVYI	VMEL	MDANL	CQVIQ	120
121	MELDHER	MSY	LLYQML	CGIK	HLHSAG	IHR	DLKPS	NIVVK	SDCTL	KILDF	GLART	AGTSF	180
181	MMPYVV	TRY	YRAPE	VILGM	GYKEN	VDLWS	VGCIM	GEMVC	HKILF	PGRDY	IDQWN	KVIEQ	240
241	LGTCP	PEFMK	KLQPT	VRTYV	ENRPKY	AGYS	FEKLF	PDVLF	PADSEH	NKLK	ASQAR	DLLSK	300
301	MLVIDA	SKRI	SVDEA	LQHPY	INVWYD	PSEA	EAPPPK	IPDK	QLDERE	HTIE	EWKELI	YKEV	360
361	MDLEER	TKNG	VIRGQ	PSPLA	QVQQ								420

blue: JNK1 sequence expressed in recombinant protein

¹[NCBI/Protein](#) accession number NP_002741.1