

Certificate of Analysis



JNK1 (non activated)

c-Jun N-terminal kinase 1

Recombinant Human Protein Kinase

HGNC Symbol: MAPK8

Synonyms: MAPK8, PRKM8, SAPK1

Product No.: 0458-0000-2

Lot: 006

Description: Human JNK1, full length, amino acids M₁-Q₃₈₄ (as in NCBI/Protein entry NP_002741.1), not activated, N-terminal HIS₆ fusion protein with a TEV cleavage site, expressed in E.coli

Product identity: JNK1 Lot 006, was confirmed as JNK1 with a JNK1 specific Western blot

Theoretical MW_{Fusion Protein}: 46,410 Da

Expression: E.coli

Purification: Immobilized Metal Affinity Chromatography

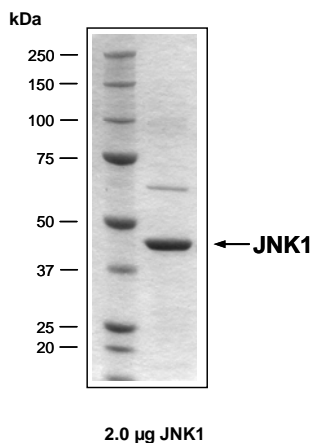
Activation: no activation

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

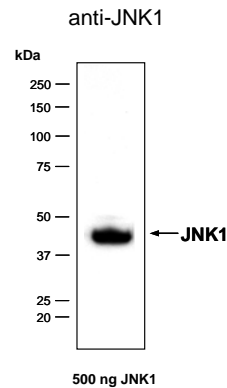
Storage temperature: -80°C
Avoid repeated freeze-thaw cycles!

Protein concentration: 0.420 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

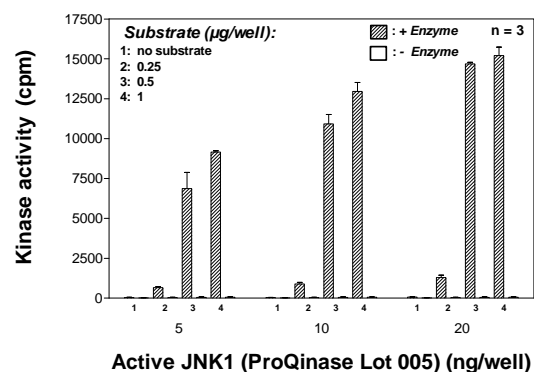
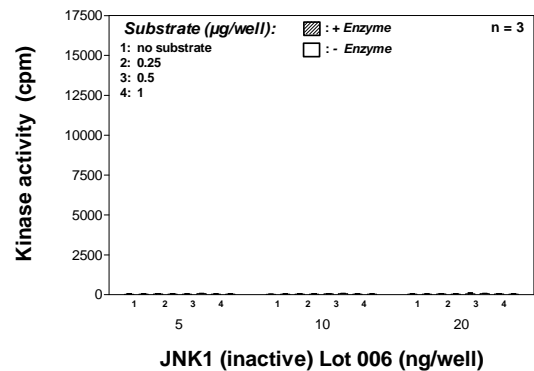
JNK1 Lot 006:
Coomassie stain



JNK1 Lot 006:
Western Blot Analysis



Kinase activity JNK1 (inactive) vs active JNK1:



Final assay concentrations:

- 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 (610,000 cpm ³³P-γ-ATP)
 - Substrate (variable): ATF2
 - Recombinant JNK1 (inactive) or active JNK1 (variable)
- Assay:** ³³PanQinase® Assay

This product is for in vitro research use only, not for use in humans or animals. ProQinase disclaims any warranty explicitly or implied that the use of the product or parts of the product is free from third party intellectual property claims unless this is explicitly stated.

CoA V1.0 JNK1_Lot006_V3.doc

JNK1 (non activated)

Product No.: 0458-0000-2

JNK1 Recombinant Fusion Protein Amino Acid Sequence														
1	M	HHHHHH	AMT	ENLYFQ	GAMS	RSKRDN	NFY	VEIGD	STFTV	LKRYQ	NLKPI	GSGAQ	GIVCA	60
61	AYDAI	LERNV	AIKKL	SRPFQ	NQTHA	KRAYR	ELVLM	KCVNH	KNIIG	LLNVF	TPQKS	SLEEFQ	120	
121	DVYIV	MELMD	ANLCQ	VIQME	LDHER	MSYLL	YQMLC	GIKHL	HSAGI	IHRDL	KPSNI	VVKSD	180	
181	CTLKI	LDFGL	ARTAG	TSFMM	TPYVV	TRYR	APEVI	LGMGY	KENVD	LWSVG	CIMGEM	VCHK	240	
241	ILFPGR	DYID	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	DEREHT	IEEW	KELIY	KEVMD	LEERT	KNGVI	RGQPS	PLAQV	QQ				420	

1-218: GST Red: HIS6-tag Green: TEV blue: JNK1

JNK1 wt ¹ amino acid sequence													
1	MSRSKR	DNNF	YSVEI	GDSTF	TVLKRY	QNLK	PIGSGA	QGIV	CAAYDA	ILERNV	NVAIKK	LSRP	60
61	FQNQTH	AKRA	YRELV	LKCV	NHKNI	IGLLN	VFTPQK	SLEE	FQDVYI	VMEL	MDANLC	QVIQ	120
121	MELDHER	MSY	LLYQML	CGIK	HLHSAG	IHR	DLKPSN	IVVK	SDCTLK	ILDF	GLARTAG	TSF	180
181	MMTPYV	VTRY	YRAPE	VILGM	GYKEN	VDLWS	VGCIMG	EMVC	HKILFP	GRDY	IDQWNK	VIEQ	240
241	LGTPCPE	FMK	KLQPT	VRTYV	ENRPKY	AGYS	FEKLPD	VLF	PADSEH	NKLN	ASQARD	LLSK	300
301	MLVIDAS	KRI	SVDEA	LQHPY	INVWYD	PSEA	EAPPPK	IPDK	QLDERE	HTEI	EWKELI	YKEV	360
361	MDLEERT	KNG	VIRGQ	PSPLA	QVQQ								420

blue: JNK1 sequence expressed in fusionprotein

¹NCBI/Protein accession number NP_002741.1