

ProQinase™ JNK2 (non activated)

mitogen-activated protein kinase 9

Recombinant Human Protein Kinase

Synonyms: MAPK9, JNK2A, JNK2ALPHA, JNK2B, JNK2BETA, JNK-55, p54a, p54aSAPK, PRKM9, SAPK, JNK2-a2

Product No.: 0459-0000-2

Lot: 004

Description: Human JNK2, full length, amino acids M₁-R₄₂₄ (as in NCBI/Protein entry NP_002743.3), N-terminal HIS₆ fusion protein with a TEV cleavage site, expressed in E.coli

Theoretical MW: 51,666 Da

Expression: E.coli

Purification: Immobilized Metal Affinity Chromatography

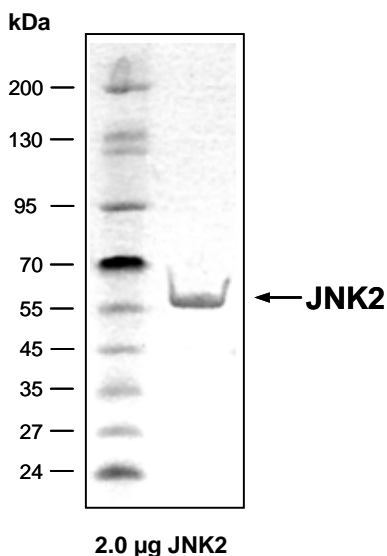
Storage buffer: 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 20% glycerol

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

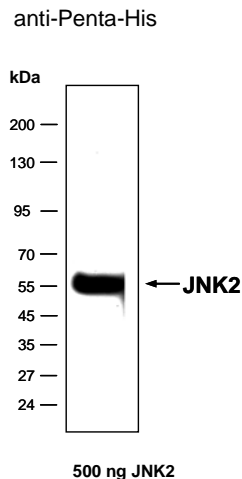
Protein concentration: 0.763 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

JNK2 Lot 004

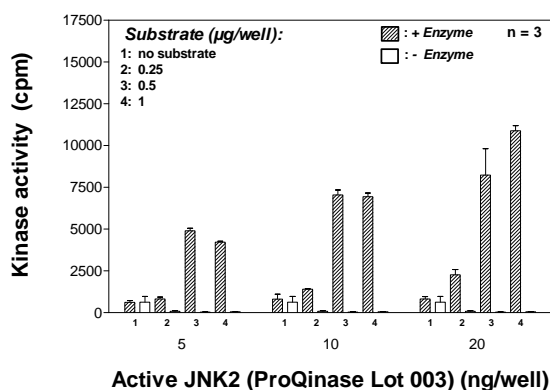
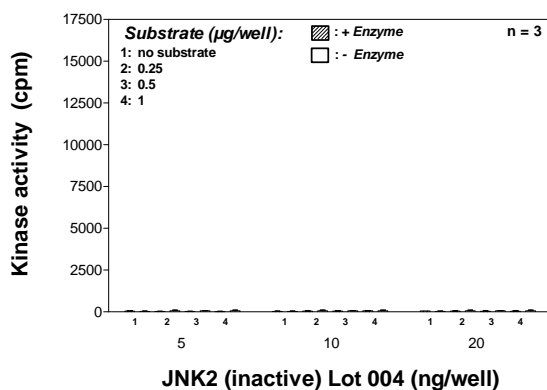
Coomassie stain:



JNK2 Lot 004: Western Blot Analysis



Kinase activity JNK2 (inactive) vs active JNK2:



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 (690,000 cpm ³³P-γ-ATP)
- Substrate (variable): ATF2
- Recombinant JNK2 (inactive) or active JNK2 (variable)

Assay: ³³PanQinase® Assay

ProQinase™ JNK2

Product No.: 0459-0000-2

JNK2 Recombinant Fusion Protein Amino Acid Sequence														
1	M	HHHHHH	AMT	ENLYFQ	GAMS	DSKCD	SQFYS	VQVAD	STFTV	LKRYQ	QLKPI	GSGAQ	GIVCA	60
61	AFD	TVLGINV	AVK	LSRPFQ	NQTHA	KRAYR	ELVLL	KCVNH	KNIIS	LLNVF	TPQKT	LEEFQ	120	
121	DVYL	VMELMD	ANLC	QVIHME	LDHER	MSYLL	YQMLC	GIKHL	HSAGI	IHRDL	KPSNI	VVKS	180	
181	CTLK	ILDFGL	ARTAC	TNFMM	TPYVV	TRYR	APEVI	LGMGY	KENVD	IWSVG	CIMGEL	VKGC	240	
241	VIFQ	TDHID	QWNK	VIEQLG	TPSAE	FMKKL	QPTVR	NYVEN	RP	RYPGI	KIFE	ELFPD	WIFPS	300
301	ESER	DKIKTS	QARDL	LSKML	VIDPD	KRISV	DEALR	HPYIT	VWYDP	AEAEA	PPPQI	YDAQL	360	
361	EERE	HAIEEW	KELIY	KEVMD	WEERS	KNGVV	KDQPS	DAAVS	SNATP	SQSSS	INDISS	MSTE	420	
421	QTLA	SDTDSS	LDAST	GPLEG	CRGK	GEFQHT	GGRY						480	

Red: HIS6-tag Green: TEV blue:JNK2 R: K265R variation

JNK2 wt ¹ ORF (¹ NCBI/Protein accession number NP_002743.3)														
1	MSD	SKCD	SQF	YSVQ	VADSTF	TVLKRY	QQLK	PIGSG	AQGIV	CAAFD	TVLGI	NVAVK	LSRP	60
61	FQN	QTHAKRA	YREL	VLLKCV	NHKNI	IISLLN	VFTPQ	KTLEE	FQDVY	LMEL	MDANL	CQVIH	120	
121	MEL	DHERMSY	LLYQ	MLCGIK	HLHS	AGIIHR	DLKPS	NIVVK	SDCTL	KILDF	GLART	ACTNF	180	
181	MMTP	YVVTRY	YRAPE	VILGM	GYKEN	VDIWS	VGCIM	GELVK	GCVIF	QGDH	IDQWN	KVIEQ	240	
241	LGTP	SAEFMK	KLQPT	VRNYV	ENRP	KYPGIK	FEELF	PDWIF	PSESE	RDKIK	TSQAR	DLLSK	300	
301	MLVI	DPDKRI	SVDEA	LHPY	ITVWY	DPAEA	EAPPP	QIYDA	QLEER	EHAIE	EWKEL	IYKEV	360	
361	MDWE	ERSKNG	VVKDQ	PSDAA	VSSNA	TSPQS	SSINDI	SSMS	TEQTL	ASDTD	SSLD	ASTGPL	420	
421	EGCR												480	

blue: sequence of JNK2 expressed in fusionprotein RED: K265 variation site

Recombinant Proteins