

JNK3 (non activated)

c-JUN N-terminal kinase 3

Recombinant Human non activated Protein Kinase

HGNC Symbol: MAPK10

Synonyms: JNK3A, p54bSAPK, PRKM10

Product No.: 0900-0000-2

Lot: 005

Description: Human JNK3, full length, amino acids M₁-R₄₂₆ (as in NCBI/Protein NP_620446.1), non activated, untagged, expressed in E.coli

Theoretical MW_{Protein}: 48,467 Da

Expression: E. coli

Purification: GST-Affinity Chromatography followed by cleavage of GST-tag

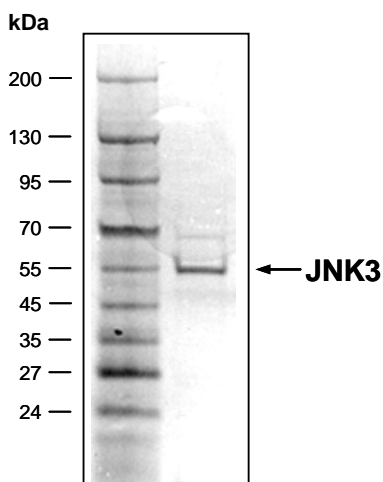
Activation: No special measures were taken to activate JNK3 Lot 005

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!

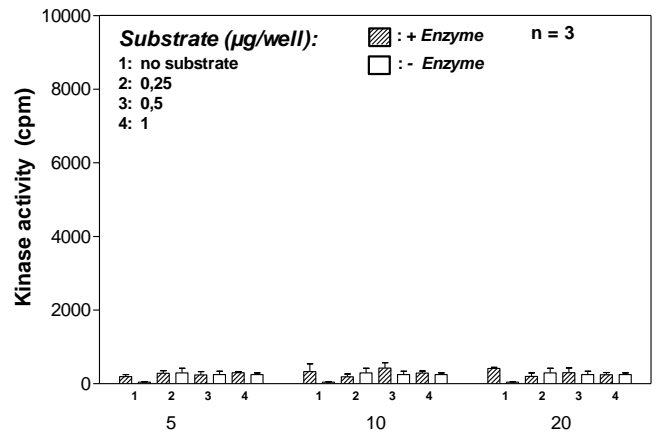
Protein concentration: 0.107 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Coomassie stain:

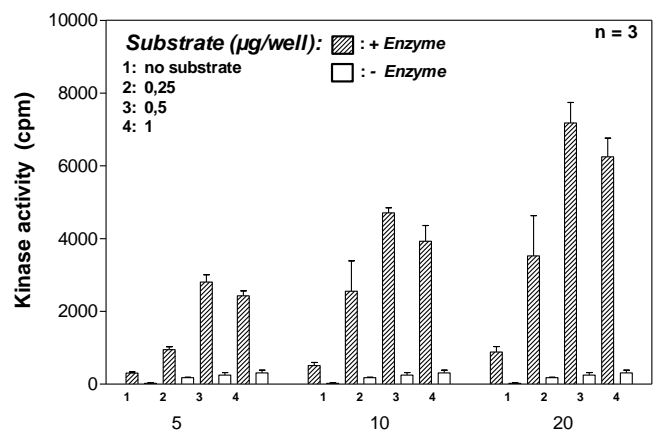


2.0 µg JNK3

Kinase activity JNK3 (non activated) vs active JNK3:



JNK3 (non activated) Lot 005 (ng/well)



Active JNK3 Lot 004 (ng/well)

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 JNK3 (non activated) or active JNK3 (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.

JNK3 (non activated)

Product No.: 0900-0000-2

JNK3 Recombinant Fusion Protein Amino Acid Sequence								
1	GPLA	MSKSKV	DNQFYSVEVG	DSTFTVLKRY	QNLKPIGSGA	QGIVCAAYDA	VLDNRNVAIKK	60
61	LSRPFQ	NQTH	AKRAYRELVL	MKCVNHKNII	SLLNVFTPOK	TLEEFQDVYL	VMELMDANLC	120
121	QVIQ	MELDHE	RMSYLLYQML	CGIKHLHSAG	IIHRDLKPSN	IVVKS DCTLK	ILDFGLARTA	180
181	GTSF	MMTPYV	VTRYRAPEV	ILGMGYKENV	DIWSVGCIMG	EMVRHKILFP	GRDYIDQWNK	240
241	VIEQLG	TPCP	EFMKKLQPTV	RNYVENRPKY	AGLTFPKLFP	DSLFPADSEH	NKLNKASQARD	300
301	LLSK	MLVIDP	AKRISVDDAL	QHPYINVWYD	PAEVEAPPPQ	IYDKQLDERE	HTIEEWKELI	360
361	YKEVMN	SEEK	TKNGVVKGQP	SPSGAAVNSS	ESLPPSSSVN	DISSMSTDQT	LASDTDSSLE	420
	ASAG	PLGCCR						420

1-4: remnant from 3C cleavage **blue: JNK3**

JNK3 wt ¹ amino acid sequence										
1	MSKSKVDNQF	YSVEVGDSTF	TVLKRYQNLK	PIGSGAQGIV	CAAYDAVLDR	NVAIKKLSRP	60			
61	FQNT	HAKRA	YRELVLKCV	NHKNIISLLN	VFTPOK	TLLEE	FQDVYLMEL	MDANLCQVIQ	120	
121	MELD	HERMSY	LLYQMLCGIK	HLHSAGIIHR	DLKPSNIVVK	SDCTLKILDF	GLARTAGTSF	180		
181	MMTPYV	VTRY	YRAPEVILGM	GYKENVDIWS	VGCIMGEMVR	HKILFPGRDY	IDQWNKVIEQ	240		
241	LGTPC	PEFMK	KLQPTVRNYV	ENRPKYAGLT	FPKLPD	SLF	PADSEHNKLNK	ASQARDLLSK	300	
301	MLVID	PAKRI	SVDDALQHPY	INVWYDPAEV	EAPPPQ	IYDK	QLDEREHTIE	EWKELIYKEV	360	
361	MNSEEK	TKNG	VVKGQPSPSG	AAVNSS	ESLP	PSSSVN	DISS	MSTDQTLASD	TDSSLEASAG	420
421	PLGCCR							480		

blue: JNK3 sequence expressed in fusionprotein

¹NCBI/Protein accession number NP_620446.1