

ProQinase™ 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

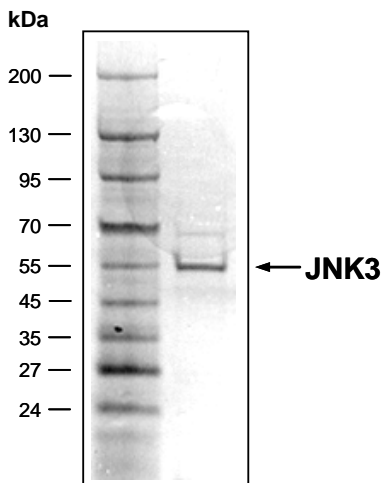
Activation: No special measures were taken to activate JNK3 Lot 005

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!

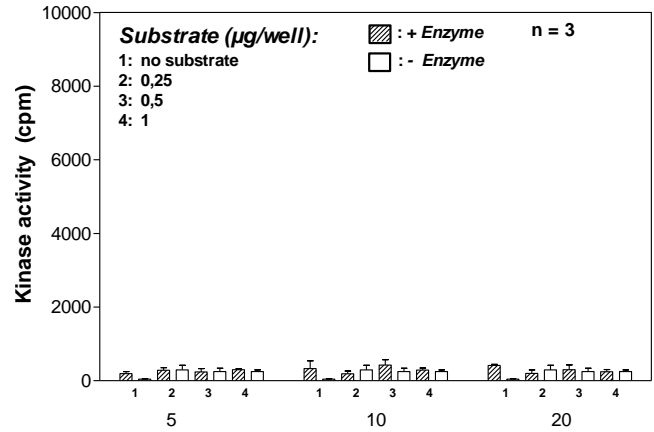
Protein concentration: 0.107 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Coomassie stain:

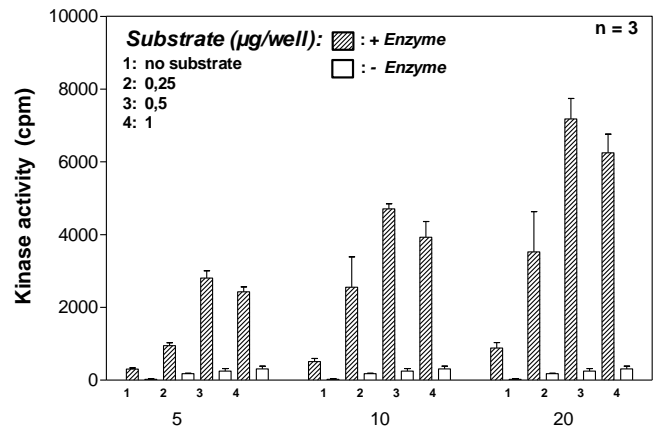


2.0 µg JNK3

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

Recombinant Proteins

ProQinase™ JNK3 (non activated)

Product No.: 0900-0000-2

JNK3 Recombinant Fusion Protein Amino Acid Sequence					
1	GPI	AMSKSKV	DNQFYSVEVG	DSTFTVLKRY QNLKPIGSGA QGIVCAAYDA VLDRNVAIKK	60
61	LSRPFQ	NQTH	AKRAYRELVL	MKCVNHKNII SLLNVFTPQK TLEEFQDVYL VMELMDANLC	120
121	QVIQ	MELDHE	RMSYLLYQML	CGIKHLHSAG I IHRDLKPSN IVVKS DCTLK ILDFGLARTA	180
181	GTSFM	MTPYV	VTRY	YRAPEV ILGMGYKENV DIWSVGCIMG EMVRHKILFP GRDYIDQWNK	240
241	VIEQL	GTPCP	EFMKKLQPTV	RNYVENRPKY AGLTFPKLFP DSLFPADSEH NKLKASQARD	300
301	LLSKM	LVIDP	AKRISVDDAL	QHPYINVWYD PAEVEAPPPQ IYDKQLDERE HTIEEWKELI	360
361	YKEVM	NSEEK	TKNGVVKGQP	SPSGAAVNSS ESLPPSSSVN DISSMSTDQT LASDTSLSLE	420
	ASAG	PLGCCR			

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

JNK3 wt ¹ amino acid sequence					
1	MSKSKV	DNQF	YSVEVGDSTF	TVLKRYQNLK PIGSGAQGIV CAAYDAVLDR NVAIKKLSRP	60
61	FQNQ	THAKRA	YRELVLKCV	NHKNIISLLN VFTPQK TLEE FQDVYLMEL MDANLCQVIQ	120
121	MELD	HERMSY	LLYQMLCGIK	HLHSAGIIHR DLKPSNIVVK SDCTLKILDF GLARTAGTSF	180
181	MTPY	VVTRY	YRAPEVILGM	GYKENVDIWS VGCIMGEMVR HKILFPGRDY IDQWNKVIEQ	240
241	LGTP	CPEFMK	KLQPTVRNYV	ENRPKYAGLT FPKLFPDSLFPADSEHNKLK ASQARDLLSK	300
301	MLVID	PAKRI	SVDDALQHPY	INVWYDPAEV EAPPPQIYDK QLDEREHTIE EWKELIYKEV	360
361	MNSEEK	TKNG	VVKGQPSPSG	AAVNSSSESLP PSSSVNDISS MSTDQTLASD TSSLEASAG	420
421	PLGCCR				480

blue: JNK3 sequence expressed in fusionprotein

¹NCBI/Protein accession number NP_620446.1

Recombinant Proteins