

ProQinase™ JUN

Jun proto-oncogene, AP-1 transcription factor subunit

Recombinant Protein Kinase Substrate

HGNC Symbol: JUN

Synonyms: AP-1, c-Jun

Product No.: 0718-0000-1

Lot: 005

Description: Human JUN, N-terminal fragment, amino acids E₅-T₈₉ (as in [NCBI/Protein](#) entry NP_002219.1), N-terminal GST-HIS₆ fusion protein with a Thrombin cleavage site, expressed in E.coli

Theoretical MW_{Fusion Protein}: 36,096 Da

Expression host: E.coli

Purification: GST-Affinity Chromatography

ATPase activity: In an ADP-Glo™ assay (Promega) with 10 μM ATP or 30 μM ATP, the ATP → ADP conversion within 30 min is approx. 1% at a concentration of 100 μg/ml substrate.
Detailed ATPase assay conditions on request

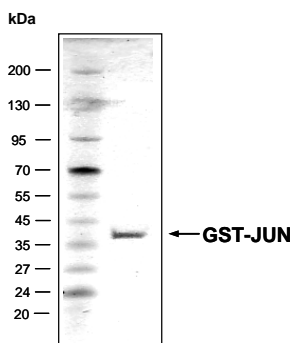
Storage buffer: 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, 20 % glycerol

Storage temperature: -80°C

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

Protein concentration: .1257 μg/μl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

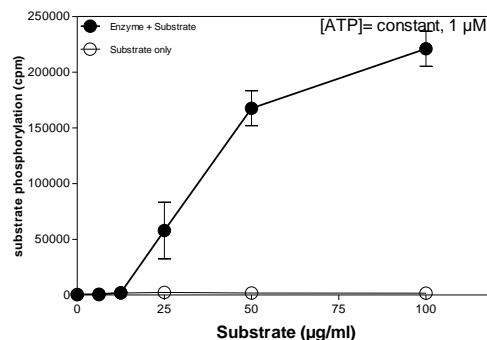
JUN Lot 005:
Coomassie stain



2.0 μg GST-JUN

Phosphorylation of JUN by the kinase JNK1

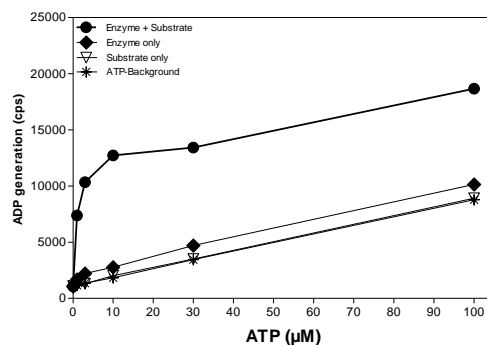
Radiometric filter binding assay



Assay conditions:

70 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}
ATP: 1 μM
Substrate (JUN): variable concentration
JNK1: 200 ng/ml
MSFC membrane (Millipore)

ADP-Glo™ assay (Promega)



70 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}
ATP: variable concentration
1% (v/v) DMSO
Substrate (JUN): 100 μg/ml
JNK1: 200 ng/ml

This product was manufactured at Reaction Biology in Freiburg, Germany, and is for in vitro research use only, not for use in humans or animals.
© European Union, 2020. Material may not be reproduced or distributed without written permission from Reaction Biology Europe GmbH

ProQinase™ JUN

Product No.: 0781-0000-1

GST-JUN Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP EML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPQID	KYLKSSKYIA	WPLQGWAQATF	GGGDHPPKSD	LVPRGSETTF	YDDALNASFL	240
241	PSESGPYGYS	NPKILKQSM T	LNLADPVGSL	KPHLRAKNSD	LLTSPDVGLL	KLASPELERL	300
301	LIQSSNGHIT	EFIVTD					360

1-218: GST **Pink**: Thrombin cleavage site **blue**: JUN fragment

This product was manufactured at Reaction Biology in Freiburg, Germany, and is for in vitro research use only, not for use in humans or animals.
© European Union, 2020. Material may not be reproduced or distributed without written permission from Reaction Biology Europe GmbH