

IKK-epsilon

inhibitor of nuclear factor kappa-B kinase epsilon subunit

Recombinant Human Active Protein Kinase

HGNC Symbol: IKBKE

Synonyms: IKK-I; IKKE; IKKI

Product No.: 0320-0000-1

Lot: 007

Description: Human IKK-epsilon, full length, amino acids M₁-V₇₁₆ (as in [NCBI/Protein](#) entry NP_054721.1), N-terminal GST-HIS₆ fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

Product identity: IKK-epsilon Lot 007, was confirmed as IKK-epsilon by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 110,059 Da

Expression host: Sf9 insect cells

Purification: GST-Affinity Chromatography

Activation: This kinase was not activated by special procedures

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 10µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 0.055 µg/µl

(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

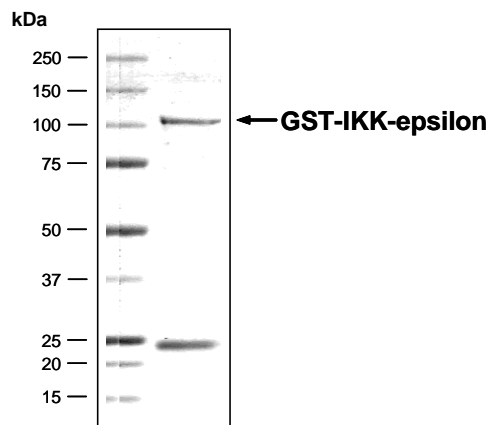
Specific kinase activity (P_i transfer): 42 pmol/µg × min

ATP-K_M: 2.3 µM

Additional assay technology:

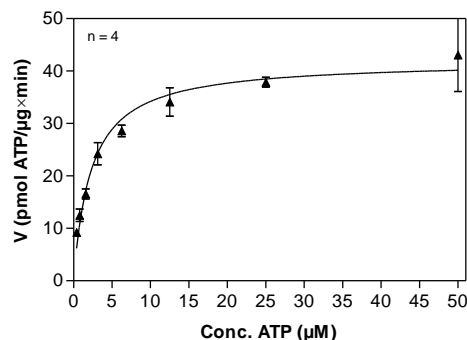
IKK-epsilon Lot 007 was also successfully tested by ProQinase for the use with the ADP-Glo™ Kinase assay from Promega. ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details

IKK-epsilon Lot 007: Coomassie stain



2.0 µg GST-IKK-epsilon

IKK-epsilon Lot 007: Determination of V_{max} and K_M value for ATP



Determination of K_M value & Specific activity:

- Assay conditions:
 - 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}
 - ATP (variable)
 - Substrate: RBER-GSK3, 100 µg/ml
 - Kinase: 1 µg/ml
- Filter binding assay
- MSFC membrane (Millipore)

IKK-epsilon

Product No.: 0320-0000-1

GST-IKK-epsilon Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPOID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHG	RRRASVAAGI	240
241	LVPRGSPGLD	GICSRNSMQS	TANYLWHTDD	LLGQGATASV	YKARNKKSQE	LVAVKVFNTT	300
301	SYLRPREVQV	REFEVLRLKN	HQNIKLVFAV	EETGGSRQKV	LVMEYCSSGS	LLSVLESPEN	360
361	AFGLPEDEFL	VVLRCSVAGM	NHLRENGIVH	RDIKPGNIMR	LVGEEGQSIY	KLTFDGAARE	420
421	LDDDEKFSVS	YGTEEYLHPD	MYERAVLRKP	QQKAFGVTVD	LWSIGVTLYH	AATGSLPFIP	480
481	FGGPRRNKEI	MYRITTEKPA	GAIAGAQRRE	NGPLEWSYTL	PITCQLSLGL	QSQLVPILAN	540
541	ILEVEQAKCW	GFDQFFAETS	DILQRVVVHV	FSLSQAVLHH	IYIHAHNTIA	IFQEAVHKQT	600
601	SVAPRHQEYL	FEGHLCVLEP	SVSAQHIAHT	TASSPLTLFS	TAIPKGLAFR	DPALDVPKRV	660
661	PKVDLQADYN	TAKGVLGAGY	QALRLARALL	DGQELMFRGL	HWVMEVLQAT	CRRTLEVART	720
721	SLLYLSSSLG	TERFSSVAGT	PEIQELKAAA	ELRSRLRTLA	EVLSRCSQNI	TETQESLSSL	780
781	NRELKSRDQ	VHEDRSIQOI	QCCLDKMFI	YKQFKKSRMR	PGLGYNEEQI	HKLDKVNFSH	840
841	LAKRLLQVFO	EECVQKYQAS	LVTHGKRMRV	VHETRNHLRL	VGCSVAACNT	EAQGVQESLS	900
901	KLLEELSHQL	LQDRAKGAQA	SPPPIAPYPS	PTRKDLLLHM	QELCEGMKLL	ASDLLDNNRI	960
961	IERLNRVPAP	PDV					1020

1-218: GST **Red**: HIS6-tag **Pink**: Thrombin cleavage site **blue**: IKK-epsilon

IKK-epsilon wt ¹ Amino Acid Sequence							
1	MQSTANYLWH	TDDLQGQAT	ASVYKARNKK	SGELVAVKVF	NTTSYLRPRE	VQVREFEVLK	60
61	KLNHQNIKLV	FAVEETGGSR	QKVLVMEYCS	SGSLLSVLES	PENAFGLPED	EFLVLRCSV	120
121	AGMNHLENG	IVHRDIKPGN	IMRLVGEEQ	SIYKLTDFGA	ARELDDDEKF	VSVYGTEEYL	180
181	HPDMYERAVL	RKPOQKAFGV	TVDLWSIGVT	LYHAATGSLP	FIPFGGPRRN	KEIMYRITTE	240
241	KPAGAIAGAQ	RRENGPLEWS	YTLPTICQLS	LGLQSQLVPI	LANILEVEQA	KCWGFDQFFA	300
301	ETSDILQRVV	VHVFSLSQAV	LHHIYIHAHN	TIAIFQEAVH	KQTSVAPRHQ	EYLFEGHLCV	360
361	LEPSVSAQHI	AHTTASSPLT	LFSTAIPKGL	AFRDPALDVP	KFVPKVDLQA	DYNTAKGVLG	420
421	AGYQALRLAR	ALLDQELMF	RGLHWVMEVL	QATCRRTLEV	ARTSLLYLSS	SLGTERFSSV	480
481	AGTPEIQELK	AAAEIERSLR	TLAEVLSRCS	QNITETQESL	SSLNRELVKS	RDQVHEDRSI	540
541	QQIQCCLDKM	NFIYKQFKKS	RMRPGLGYNE	EQIHKLDKVN	FSLAKRLLQ	VFQECCVQKY	600
601	QASLVTHGKR	MRVVHETRNH	LRLVGCSVAA	CNTEAQGVQE	SLSKLEELS	HQLLQDRAG	660
661	AQASPPPIAP	YSPTRKDLL	LHMQELCEGM	KLLASDLLDN	NRIIERLNRV	PAPPDV	720

blue: IKK-epsilon sequence expressed in recombinant protein

¹[NCBI/Protein](#) accession number NP_054721.1

This product was manufactured at ProQinase in Freiburg, Germany, and 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.