

ProQinase™ AKT1 aa106-480

AKT serine/threonine kinase 1

Recombinant Human Active Protein Kinase

HGNC Symbol: AKT1

Synonyms: PKB, RAC, PRKBA, RAC-ALPHA

Product No.: 0132-0000-2

Lot: 007

Description: Human AKT1, C-terminal fragment, amino acids V₁₀₆-A₄₈₀ (as in [NCBI/Protein](#) entry NP_005154.1), N-terminal GST-HIS₆, C-terminal HIS₆ fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

Product identity: AKT1 Lot 007, was confirmed as AKT1 by Western Blotting using a specific anti AKT1 antibody

Theoretical MW_{Fusion Protein}: 77,832 Da

Expression host: Sf9 insect cells

Purification: GST-Affinity Chromatography

Activation: With PDK1

Storage buffer: 50 mM TRIS-HCl pH 8.0, 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: 0.434 µg/µl

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

Biochemical Parameters:

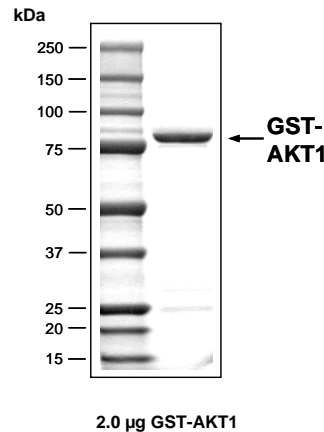
Specific kinase activity (P_i transfer): 328 pmol/µg × min

ATP-K_M: 41 µM

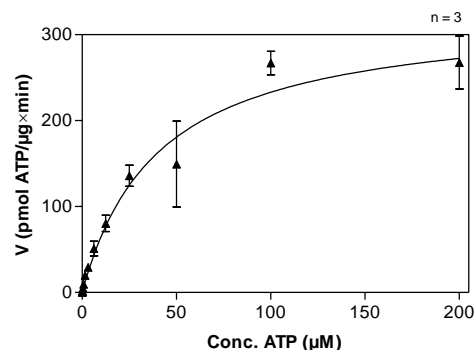
Additional assay technology:

AKT1 aa106-480 Lot 007 was also successfully tested by Reaction Biology 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

AKT1 aa106-480 Lot 007:
Coomassie stain



AKT1 aa106-480 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: R₁₁-GSK3(14-27), 100 µg/ml
 - AKT1: 2 µg/ml
- Filter binding assay
- MSPH membrane (Millipore)

ProQinase™ AKT1 aa106-480

Product No.: 0132-0000-2

GST-AKT1 aa106-480 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRLL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPOID	KYLKSSKYIA	WPLQGQWQATF	GGGDHPPKSD	PMGHHHHHG	RRRASVAAGI	240
241	LVPRGSPGLD	GIYARGIQAS	MGARGRLQGT	PVADGLKKQE	EEEMDFRSGS	PSDNSGAEEM	300
301	EVSLAKPKHR	VTMNEFEYLK	LLGKGTFGKV	ILVKEKATGR	YYAMKILKKE	VIVAKDEVVAH	360
361	TLTENRVLQN	SRHPFLTALK	YSFQTHDRLC	FVMEYANGGE	LFFHLSRERV	FSEDRARFYG	420
421	AEIVSALDYL	HSEKNVVYRD	LKLENLMLDK	DGHIKITDFG	LCKEGIKDGA	TMKTFCGTPE	480
481	YLAPEVLEDN	DYGRAVDWWG	LGVVMYEMMC	GRLPFYNQDH	EKLFELILME	EIRFPRTLGP	540
541	EAKSLLSGLL	KKDPKQRLGG	GSEDAKEIMQ	HRFFAGIVWQ	HVYEKLSPP	FKPQVTSETD	600
601	TRYFDEEFTA	QMITITPPDQ	DDSMCEVDSE	RRPHFPQFSY	SASSTAKGEL	RGHPFEGKPI	660
661	PNLLGLDST	RTGHHHHHH					720

1-218: GST **Red:** HIS6-tag **Pink:** Thrombin cleavage site **blue:** AKT1 fragment

AKT1 wt ¹ Amino Acid Sequence							
1	MSDVAIVKEG	WLHKRGEYIK	TWRPRYFLLK	NDGTFIGYKE	RPQDVDQREA	PLNNFSVAQC	60
61	QLMKTERPRP	NTFIIIRCLQW	TTVIERTFHV	ETPEEREWT	TAIQTVADGL	KKQEEEEEMDF	120
121	RSGSPSDNSG	AEEMEVS LAK	PKHRVTMNEF	EYLKLLGKGT	FGKVILVKEK	ATGRYYAMKI	180
181	LKKEVIVAKD	EVAHTLTENR	VLQNSRHPFL	TALKYSFQTH	DRLCFVMEYA	NGGELFFHLS	240
241	RERVFSEDRA	RFYGAETVSA	LDYLHSEKNV	VYRDLKLENL	MLDKDGHIKI	TDFGLCKEGI	300
301	KDGATMKTFC	GTPEYLAPEV	LEDNDYGRAV	DWWGLGVVY	EMMCGRLPFY	NQDHEKLFEL	360
361	IIMEEIRFPR	TLGPEAKSL	SGLLKKDPKQ	RLGGGSEDAK	EIMQHRFFAG	IVWQHVEYK	420
421	LSPPFKPQVT	SETDTRYFDE	EFTAQMITIT	PPDQDSMEC	VDSERRPHFP	QFSYSASSTA	480

blue: AKT1 sequence expressed in recombinant protein

¹[NCBI/Protein](https://www.ncbi.nlm.nih.gov/Protein) accession number NP_005154.1