

## AKT2 aa1-481

AKT serine/threonine kinase 2

Recombinant Human Active Protein Kinase

HGNC Symbol: AKT2

**Synonyms:** PKBB, PKBBETA, PKB beta, PRKBB, RAC-BETA, RAC-PK-beta

**Product No.:** 1577-0000-1

**Lot:** 003

**Description:** Human AKT2, full length, amino acids M<sub>1</sub>-E<sub>481</sub> (as in [NCBI/Protein](#) entry NP\_001617.1), N-terminal GST-HIS<sub>6</sub> fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

**Product identity:** AKT2 Lot 002, was confirmed as AKT2 by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW**<sub>Fusion Protein</sub>: 84,273 Da

**Expression host:** Sf9 insect cells

**Purification:** GST-Affinity Chromatography

**Activation:** With PDK1

**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:** 0.471 µg/µl

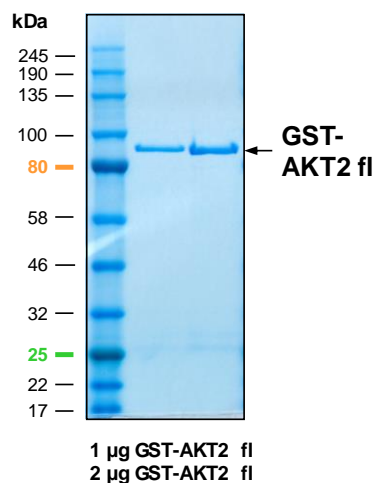
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

**Biochemical Parameters:**

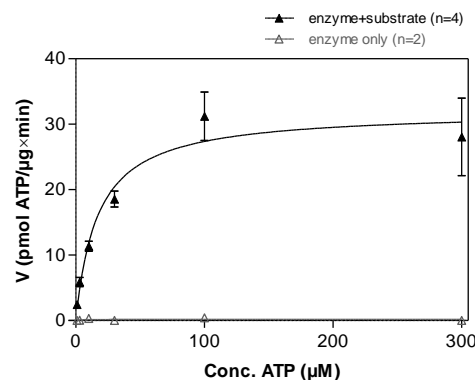
Specific kinase activity (P<sub>i</sub> transfer): 32 pmol/µg × min

ATP-K<sub>M</sub>: 17 µM

### AKT2 aa1-481 Lot003: Coomassie stain



### AKT2 aa1-481 Lot003: Determination of V<sub>max</sub> and K<sub>M</sub> value for ATP



### Determination of K<sub>M</sub> value & Specific activity:

- Assay conditions:
  - 60 mM HEPES-NaOH, pH 7.5
  - 3 mM MgCl<sub>2</sub>
  - 3 mM MnCl<sub>2</sub>
  - 3 µM Na-orthovanadate
  - 1.2 mM DTT
  - 50 µg/ml PEG<sub>20,000</sub>
  - ATP (variable)
  - Substrate: GSK3-derived peptide 100 µg/ml
  - Kinase: 1 µg/ml
- Filter binding assay
  - MSPH membrane (Millipore)

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.

## AKT2 aa1-481

Product No.: 1577-0000-1

GST-AKT2 aa1-481 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEML	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMG <b>HHHHHG</b>	RDS <b>LEVLFQG</b>	240
241	<b>PLAMVMNEVS</b>	<b>VIKEGWLHKR</b>	<b>GEYIKTWRPR</b>	<b>YFLLKSDGSF</b>	<b>IGYKERPEAP</b>	<b>DQTLPLNNEF</b>	300
301	<b>SVAECQLMKT</b>	<b>ERPRPNTFVI</b>	<b>RCLQWTTVIE</b>	<b>RTFHVDSPE</b>	<b>REEWMRAIQM</b>	<b>VANSLKQRAP</b>	360
361	<b>GEDPMDYKCG</b>	<b>SPSDSSTTEE</b>	<b>MEVAVSKARA</b>	<b>KVTMNDFDYL</b>	<b>KLLGKGTFGK</b>	<b>VILVREKATG</b>	420
421	<b>RYYAMKILRK</b>	<b>EVIIAKDEVA</b>	<b>HTVTESRVLQ</b>	<b>NTRHPFLTAL</b>	<b>KYAFQTHDRL</b>	<b>CFVMEYANGG</b>	480
481	<b>ELFFHLSRER</b>	<b>VFTEERARFY</b>	<b>GAEIVSALEY</b>	<b>LHSRDVVYRD</b>	<b>IKLENLMLDK</b>	<b>DGHIKITDFG</b>	540
541	<b>LCKEGISDGA</b>	<b>TMKTFCGTPE</b>	<b>YLAPEVLEDN</b>	<b>DYGRAVDWWG</b>	<b>LGVVMYEMMC</b>	<b>GRLPFYNQDH</b>	600
601	<b>ERLFELILME</b>	<b>EIRFPRTLSP</b>	<b>EAKSLLAGLL</b>	<b>KKDKPKQLGG</b>	<b>GPSDAKEVME</b>	<b>HRFFLSINWQ</b>	660
661	<b>DVVQKLLPP</b>	<b>FKPQVTSEVD</b>	<b>TRYFDDEFTA</b>	<b>QSITITPPDR</b>	<b>YDSLGLLELD</b>	<b>QRTHFPPQFSY</b>	720
721	<b>SASIRE</b>						780

1-218: GST **Red**: HIS6-tag **Green**: 3C cleavage site **blue**: AKT2

AKT2 wt <sup>1</sup> Amino Acid Sequence							
1	<b>MNEVSVIKEG</b>	<b>WLHKRGEYIK</b>	<b>TWRPRYFLLK</b>	<b>SDGSFIGYKE</b>	<b>RPEAPDQTLP</b>	<b>PLNNFSVAEC</b>	60
61	<b>QLMKTERPRP</b>	<b>NTFVIRCLQW</b>	<b>TTVIERTFHV</b>	<b>DSPDEREEM</b>	<b>RAIQMVANSL</b>	<b>KQAPGEDPM</b>	120
121	<b>DYKCGSPSDS</b>	<b>STTEEMEVAV</b>	<b>SKARAKVTMN</b>	<b>DFDYLLKLGK</b>	<b>GTFGKVLVLR</b>	<b>EKATGRYYAM</b>	180
181	<b>KILRKEVIAA</b>	<b>KDEVAHTVTE</b>	<b>SRVLQNRHP</b>	<b>FLTALKYAFQ</b>	<b>THDRLCFVME</b>	<b>YANGGELFFH</b>	240
241	<b>LSRERVFTTE</b>	<b>RARFYGAEIV</b>	<b>SALEYLHSRD</b>	<b>VVYRDIKLEN</b>	<b>LMLDKDGHK</b>	<b>ITDFGLCKEG</b>	300
301	<b>ISDGATMKT</b>	<b>CGTPEYLAP</b>	<b>VLEDNDYGRA</b>	<b>VDWWGLGVVM</b>	<b>YEMMCGRLPF</b>	<b>YNQDHERLFE</b>	360
361	<b>LILMEEIRFP</b>	<b>RTLSPEAKSL</b>	<b>LAGLLKKDPK</b>	<b>QRLGGGSPDA</b>	<b>KEVMEHRFFL</b>	<b>SINWQDVVQK</b>	420
421	<b>KLLPPFKPQV</b>	<b>TSEVDTRYFD</b>	<b>DEFTAQSITI</b>	<b>TPPDYDSL</b>	<b>LLELDQRTHE</b>	<b>PQFSYSASIR</b>	480
481	<b>E</b>						540

**blue**: AKT2 sequence expressed in recombinant protein

<sup>1</sup>[NCBI/Protein](#) accession number NP\_001617.1

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.