

## ProQinase™ HIPK3

homeodomain interacting protein kinase 3

Recombinant Human Active Protein Kinase

HGNC Symbol: HIPK3

Synonyms: ANPK, DYRK6, FIST, FIST3, PKY, YAK1

Product No.: 0936-0000-1

Lot: 004

**Description:** Human HIPK3, internal fragment, amino acids T<sub>163</sub>-N<sub>562</sub> (as in NCBI/Protein entry NP\_005725.3), activated, untagged, expressed in Sf9 insect cells

**Product identity:** HIPK3 Lot 004, was confirmed as HIPK3 by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW<sub>Protein</sub>:** 46,163 Da

**Expression:** Baculovirus infected Sf9 cells

**Purification:** GST-Affinity Chromatography, followed by cleavage of GST-tag

**Activation:** In vitro autoactivation

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

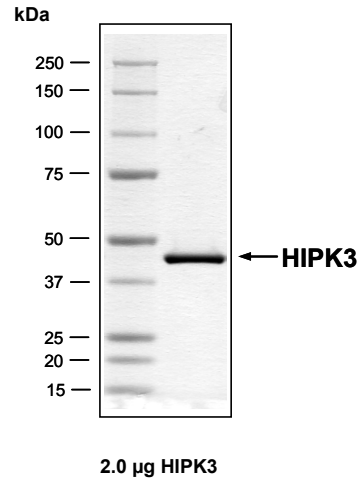
**Storage temperature:** -80°C  
Avoid repeated freeze-thaw cycles!

**Protein concentration:** 0.094 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

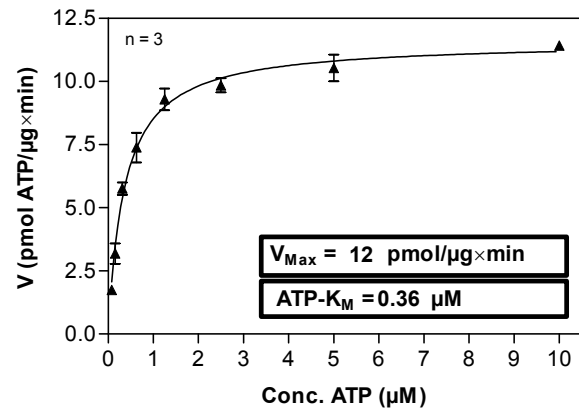
**Biochemical Parameters:**

Specific activity: 12 pmol/µg×min  
ATP-K<sub>M</sub>: 0.4 µM

**HIPK3 Lot 004:  
Coomassie stain**



**HIPK3 Lot 004:  
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: RBER-IRStide, 100 µg / ml
  - HIPK3: 1.0 µg / ml
- Filter binding assay
  - MSFC membrane (Millipore)

Recombinant Proteins

## ProQinase™ HIPK3

Product No.: 0936-0000-1

HIPK3 Recombinant Fusion Protein Amino Acid Sequence							
1	GPLAMGTVVT	ATTGSKQNC	TGEGDYQLVQ	HEVLCSMKNT	YEVLDLGRG	TFGQVVKCWK	60
61	RGTNEIVAIAK	ILKNHPSYAR	QGQIEVSILA	RLSTENADEY	NFVRAYECFQ	HRNHTCLVFE	120
121	MLEQONLYDFL	KQNKFSPLPL	KVIRPILQQV	ATALKKLSL	GLIHADLKPE	NIMLVDPVRQ	180
181	PYRVKVIDFG	SASHVSKTVC	STYLQSRYYR	APEIILGLPF	CEAIDMWSLG	CVIAELFLGW	240
241	PLYPGALEYD	QIRYISQTQ	LPGEQLLNVG	TKSTRFFCKE	TDMSHSGWRL	KTLEEHEAET	300
301	GMKSKEARKY	IFNSLDDVAH	VNTVMDLEGS	DLAEKADRR	EFVSLKML	LIDADLRITP	360
361	AETLNHPFVN	MKHLDFPHS	NHVKSCFHIM	DICKSHLNSC	DTNNHN		420

1-6: legacy from 3C cleavage    blue: HIPK3 fragment

HIPK3 wt <sup>1</sup> amino acid sequence							
1	MASQVLVYPP	YVYQTQSSAF	CSVKCLKVEP	SSCVFQERNY	PRTYVNGRNF	GNSHPPTKGS	60
61	AFQTKIPFNR	PRGHNFSLQT	SAVVLKNTAG	ATKVIAAQAQ	QAHVQAPQIG	AWRNRLLHFLE	120
121	GPQRCGLKRR	SEELDNHSSA	MQIVDELSIL	PAMLQTNMGN	PVTVVTATTG	SKQNCCTGEG	180
181	DYQLVQHEVL	CSMKNTYEVL	DFLGRGTFGQ	VVKCWKRGTN	EIVAIAKILKN	HPSYARQGQI	240
241	EVSILARLST	ENADEYNFVR	AYECFQHRNH	TCLVFEMLEQ	NLYDFLKQNK	FSPLPLKVIR	300
301	PILQQVATAL	KKLKSGLLIH	ADLKPENIML	VDPVRQPYRV	KVIDFGSASH	VSKTVCSTYL	360
361	QSRYYRAPEI	ILGLPFCEAI	DMWSLGCVIA	ELFLGWPLYP	GALEYDQIRY	ISQTQGLPGE	420
421	QLLNVGTKST	RFFCKETDMS	HSGWRKLTLE	EHEAETGMKS	KEARKYIFNS	LDDVAHVNTV	480
481	MDLEGSLLA	EKADRREFVS	LLKKMLLIDA	DLRITPAETL	NHPFVNMKHL	LDFPHSNHVK	540
541	SCFHIMDICK	SHLNSCDTNN	HNKTSLLRPV	ASSSTATLTA	NFTKIGTLRS	QALTTSAHSV	600
600	VHHGIPLQAG	TAQFGCGDAF	QQTLLIICPPA	IQGIPTHGK	PTSYSIRVDN	TVPLVTQAPA	660
661	VQPLQIRPGV	LSQTSWGRG	QMLVPAWQQV	TPLAPATTTL	TSESVAGSHR	LGDWGKMISC	720
721	SNHYNSVMPQ	PLLTNQITLS	APQPVSVGIA	HVVWPQPATT	KKNKQCQNRG	ILVKLMEWEP	780
781	GREEINAFSW	SNSLQNTNIP	HSAFISPKII	NGKDVVEVSC	IETQDNQNSE	GEARNCCETS	840
841	IRQDSDSSVS	DKQRQTIIIA	DSPSPAVSVI	TISSDTDEEE	TSQRHSLREC	KGSLDCEACQ	900
901	STLNIDRMCS	LSSPDSTLST	SSSGQSSPSP	CKRPNSMSDE	EQESSCDTVD	GSPTSDSSGH	960
961	DSPFAESTFV	EDTHENTELV	SSADTETKPA	VCSVVVPPVE	LENGLNADEH	MANTDSICQP	1020
1021	LIKGRSAPGR	LNQPSAVGTR	QQKLTSAFQQ	QHLNFSQVQH	FGSGHQEWNG	NFGHRRQQAY	1080
1081	IPTSVTSNPF	TLSHGSPNHT	AVHAHLAGNT	HLGGQPTLLP	YPSSATLSSA	APVAHLLASP	1140
1141	CTSRPMLQHP	TYNISHPSGI	VHQVPVGLNP	RLLPSPTIHQ	TQYKPIFPPH	SYIAASPAYT	1200
1201	GFPLSPTKLS	QYPYM					1260

blue: HIPK3 sequence expressed in fusionprotein

<sup>1</sup>NCBI/Protein accession number NP\_005725.3

Recombinant Proteins