

## ProQinase™ ULK3

unc-51 like kinase 3

### Recombinant Human Active Protein Kinase

HGNC Symbol: ULK3

Synonyms: DKFZP434C131; FLJ90566

Product No.: 1939-0000-1

Lot: 002

**Description:** Human ULK3, full length, amino acids M<sub>1</sub>-Q<sub>472</sub> (as in [NCBI/Protein](#) entry NP\_001092906.3), N-terminal GST-HIS<sub>6</sub> fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

**Product identity:** ULK3 Lot 002, was confirmed as ULK3 by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW**<sub>Fusion Protein</sub>: 81534 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

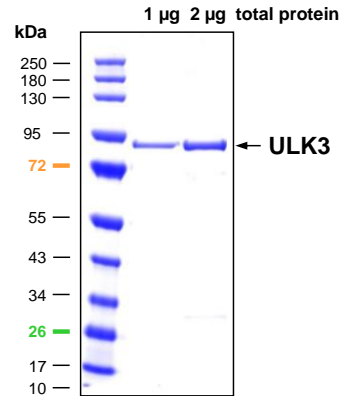
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

**Protein concentration:** 0.484 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

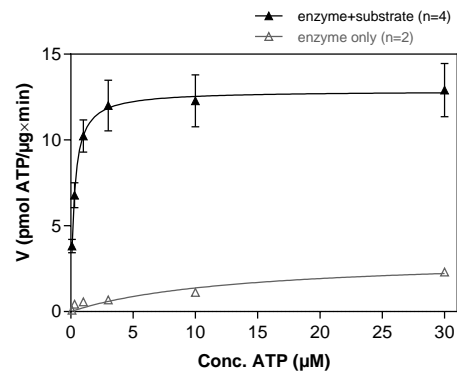
**Biochemical Parameters:**

Specific kinase activity (P<sub>i</sub> transfer): 13 pmol/µg\*min  
ATP-K<sub>M</sub>: 0.25 µM

ULK3 Lot 002:  
Coomassie stain



ULK3 Lot 002:  
Determination of V<sub>max</sub> and K<sub>M</sub> value for ATP



- 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: Casein 100 µg/ml
  - Kinase: 1 µg/ml
- Filter binding assay  
MSFC membrane (Millipore)

Recombinant Proteins

Sequence information

GST-ULK3 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRLL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEML	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPIQID	KYLKSSKYIA	WPLQGQWQATF	GGGDHPPKSD	PMGHHHHHGG	RDSLEVLFGG	240
241	MAGPGWGPP	LDGFFILTER	LGSPTYATVY	KAYAKDTRE	VVAIKCVAKK	SLNKASVENL	300
301	LTEIEILKGI	RHPHIVQLKD	FQWSDNIYL	IMEFCAGGDL	SRFIHTRRIL	PEKVARVFMQ	360
361	QLASALQFLH	ERNISHLDLK	QNILSSLE	KPHLKLADFG	FAQHMSPWDE	KHVLGRSPLY	420
421	MAPEMVCQRQ	YDARVDLWSM	GVILYEALFG	QPPFASRSFS	ELEEKIRSNR	VIELPLRPLL	480
481	SRDCRDLQRL	LLERDPSRRI	SFQDFFAHPW	VDLEHMPSGE	SLGRATALVV	QAVKKDQEGD	540
541	SAAALSLYCK	ALDFFVPALH	YEVDAQRKEA	IKAKVGQYVS	RAEELKAIVS	SSNQALLRQG	600
601	TSARDLLREM	ARDKPRLLAA	LEVASAAMAK	EEAAGGEQDA	LDLYQHSLGE	LLLLLAAEPP	660
661	GRRRELLHTE	VQNLARAAY	LKEQVKMRES	RWEADTLDE	GLSESVRSSC	TLQ	720

1-218: GST **Red**: HIS6-tag **Green**: 3C cleavage site **blue**: ULK3

ULK wt <sup>1</sup> Amino Acid Sequence							
1	MAGPGWGPPR	LDGFFILTERL	GSGTYATVYK	AYAKKDTREV	VAIKCVAKKS	LNKASVENLL	60
61	TEIEILKGI	RHPHIVQLKDF	QWSDNIYLI	MEFCAGGDLS	RFIHTTRILP	EKVARVFMQQ	120
121	LASALQFLHE	RNISHLDLKP	QNILSSLEK	PHLKLADFGF	AQHMSPWDEK	HVLGRSPLYM	180
181	APEMVCQRQY	DARVDLWSMG	VILYEALFGQ	PPFASRSFSE	LEEKIRSNRV	IELPLRPLLS	240
241	RDCRDLQRL	LERDPSRRIS	FQDFFAHPWV	DLEHMPSGES	LGRATALVVQ	AVKKDQEGDS	300
301	AAALSLYCKA	LDFFVPALHY	EVDAQRKEAI	KAKVGQYVSR	AEELKAIVSS	SNQALLRQGT	360
361	SARDLLREMA	RDKPRLLAAL	EVASAAMAKE	EAAGGEQDAL	DLYQHSLGEL	LLLLLAAEPPG	420
421	RRRELLHTEV	QNLARAAYL	KEQVKMRESR	WEADTLDEK	LSESVRSSCT	LQ	480

**blue**: ULK3 sequence expressed in recombinant protein

<sup>1</sup>[NCBI/Protein](https://www.ncbi.nlm.nih.gov/Protein) accession number NP\_001092906.3