

## ITK

IL2 inducible T cell kinase

Recombinant Human Active Protein Kinase

HGNC Symbol: ITK

Synonyms: EMT; LYK; PSCTK2

Product No.: 0359-0000-1

Lot: 001

**Description:** Human ITK, full length, amino acids M1-L<sub>630</sub> (as in [NCBI/Protein](#) entry NP\_005537.3), N-terminal GST-HIS<sub>6</sub> fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

**Product identity:** ITK Lot 001, was confirmed as ITK by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW**<sub>Fusion Protein</sub>: 101,727 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

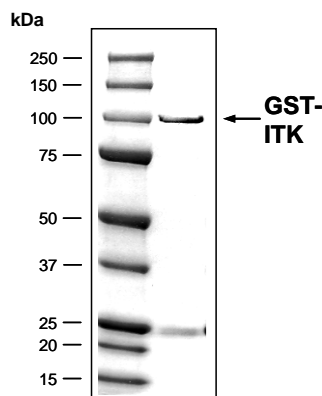
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

**Protein concentration:** 0.073 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

### Biochemical Parameters:

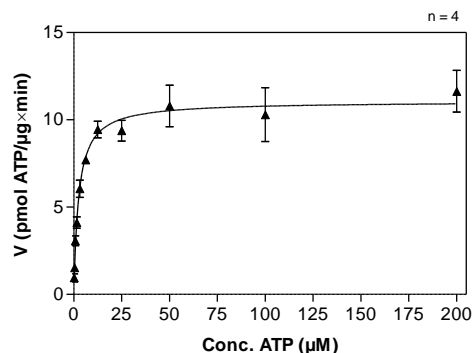
Specific kinase activity (P<sub>i</sub> transfer): 11 pmol/µg × min  
ATP-K<sub>M</sub>: 2.5 µM

### ITK Lot 001: Coomassie stain



2.0 µg GST-ITK

### ITK Lot 001: 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: Poly(Glu:Tyr)<sub>4:1</sub>, 20 µg/ml  
Kinase: 4 µg/ml
- Filter binding assay  
MSFC membrane (Millipore)

# ITK

Product No.: 0359-0000-1

GST-ITK Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSM	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHG	RRRASVAAGI	240
241	LVPRG	GIYARGIQAS	MNNFILLEEQ	LIKKSQQR	TSPSNFKVRF	FVLTKASLAY	300
301	FEDRHGKRT	LKGSIELSRI	KCVEIVKSDI	SIPCHYKYPF	QVVDNYLLY	VFAPDRESRQ	360
361	RWVLALKEET	RNNNSLVPKY	HPNFWMGKW	RCCSQLEKLA	TGCAQYDPTK	NASKKPLPPT	420
421	PEDNRRPLWE	PEETVVIALY	DYQTNDPQEL	ALRRNEEYCL	LDSSIEHWWR	VQDRNGHEGY	480
481	VPSSYLVEKS	PNNLETYEWY	NKSISRDKAE	KLLLDTGKEG	AFMVRDSRTA	GTYTTSVFTK	540
541	AVVSENNPCI	KHYHIKETND	NPKRYVVAEK	YVFDSIPLLI	NYHQHNGGGL	VTRLRYPVCF	600
601	GRQKAPVTAG	LRYGKVIDP	SELTFFVQEI	SGQFGLVHLG	YWLNKDKVAI	KTIREGAMSE	660
661	EDFIEEAEM	MKLSHPKLVQ	LYGVCLEQAP	ICLVFEMEH	GCLSDYLRTQ	RGLFAAETLL	720
721	GMCLDVCEGM	AYLEEACVIH	RDLAARNCLV	GENQVIKUSD	FGMTRFVLD	QYTSSTGTFK	780
781	PVKWASPEVF	SFSRYSSKSD	VWSFGVLMWE	VFSEGIKPYE	NRSNSEVVED	ISTGFRLYKP	840
841	RLASTHVYQI	MNHCWKERPE	DRPAFSRLLR	QLAEIAESGL			900

1-218: GST Red: HIS6-tag Pink: Thrombin cleavage site blue: ITK

ITK wt <sup>1</sup> Amino Acid Sequence							
1	MNNFILLEEQ	LIKKSQQR	TSPSNFKVRF	FVLTKASLAY	FEDRHGKRT	LKGSIELSRI	60
61	KCVEIVKSDI	SIPCHYKYPF	QVVDNYLLY	VFAPDRESRQ	RWVLALKEET	RNNNSLVPKY	120
121	HPNFWMGKW	RCCSQLEKLA	TGCAQYDPTK	NASKKPLPPT	PEDNRRPLWE	PEETVVIALY	180
181	DYQTNDPQEL	ALRRNEEYCL	LDSSIEHWWR	VQDRNGHEGY	VPSSYLVEKS	PNNLETYEWY	240
241	NKSISRDKAE	KLLLDTGKEG	AFMVRDSRTA	GTYTTSVFTK	AVVSENNPCI	KHYHIKETND	300
301	NPKRYVVAEK	YVFDSIPLLI	NYHQHNGGGL	VTRLRYPVCF	GRQKAPVTAG	LRYGKVIDP	360
361	SELTFFVQEI	SGQFGLVHLG	YWLNKDKVAI	KTIREGAMSE	EDFIEEAEM	MKLSHPKLVQ	420
421	LYGVCLEQAP	ICLVFEMEH	GCLSDYLRTQ	RGLFAAETLL	GMCLDVCEGM	AYLEEACVIH	480
481	RDLAARNCLV	GENQVIKUSD	FGMTRFVLD	QYTSSTGTFK	PVKWASPEVF	SFSRYSSKSD	540
541	VWSFGVLMWE	VFSEGIKPYE	NRSNSEVVED	ISTGFRLYKP	RLASTHVYQI	MNHCWKERPE	600
601	DRPAFSRLLR	QLAEIAESGL					660

blue: ITK sequence expressed in recombinant protein

<sup>1</sup>[NCBI/Protein](#) accession number NP\_005537.3

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.