

IRAK4 (GST-HIS-tag)

interleukin-1 receptor associated kinase 4

Recombinant Human Active Protein Kinase

HGNC Symbol: IRAK4

Synonyms: REN64; NY-REN-64; IPD1

Product No.: 0268-0000-1

Lot: 005

Description: Human IRAK4, C-terminal fragment, amino acids A₁₀₄-S₄₆₀ (as in [NCBI/Protein](#) entry NP_057207.2), N-terminally fused to GST-HIS₆ and a Thrombin cleavage site, expressed in Sf9 insect cells

Product identity: IRAK4 Lot 005, was confirmed as IRAK4 by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 69,898 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, 4 mM reduced glutathione, 20 % glycerol

Storage temperature: -80°C

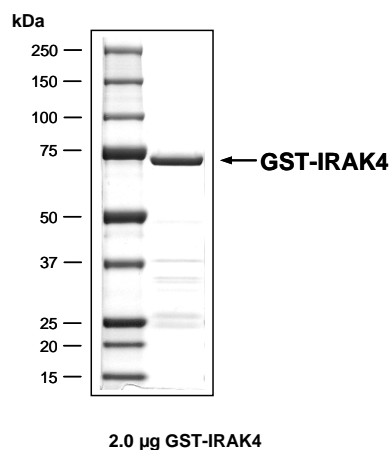
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 0.213 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

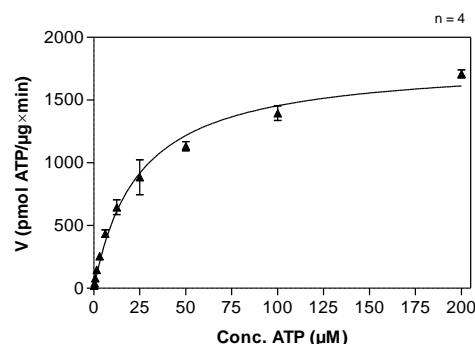
Biochemical Parameters:

Specific kinase activity (P_i transfer): 1809 pmol/µg × min
ATP-K_M: 24 µM

IRAK4 Lot 005: Coomassie stain



IRAK4 Lot 005: 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: Histon H2B 100 µg/ml
 - Kinase: 0.4 µg/ml
- Filter binding assay
MSFC 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.

IRAK4 (GST-HIS-tag)

Product No.: 0268-0000-1

GST-IRAK4 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP EML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI PQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHG	RRRASVAAGI	240
241	LVPRG SGLD	GICSIEEFTM	AVPKTANTLP	SKEAITVQOK	QMPFCDKDRT	LMPVQNL EQ	300
301	SYMPDSSSP	ENKSLEVSDT	RFHSFSFYEL	KNVTNNFDER	PISVGGNKMG	EGGFGVVYKG	360
361	YVNNTTVAVK	KLAAMVDITT	EELKQQFDQE	IKVMAKQHE	NLVELLGFSS	DGDDLCLVYV	420
421	YMPNGSLLDR	LSCLDGTPL	SWHMRCKIAQ	GAANGINFLH	ENHHIHRDIK	SANILLDEAF	480
481	TAKISDFGLA	RASEKFAQTV	MTSRIVGTTA	YMAPEALRGE	ITPKSDIYSF	GVVLEIITG	540
541	LPAVDEHREP	QLLLDIKEEI	EDEEKTIEDY	IDKKNMDADS	TSVEAMYSVA	SQCLHEKKNK	600
601	RPDIKKVQQL	LQEMTAS					660

1-218: GST **Red:** HIS6-tag **Pink:** Thrombin cleavage site **blue:** IRAK4 fragment

IRAK4 wt ¹ Amino Acid Sequence							
1	MNKPIPTSTY	VRCLNVGLIR	KLSDFIDPQE	GWKKLAVAIAK	KPSGDDRYNQ	FHIRRFEALL	60
61	QTGKSPTSEL	LFDWGTTNCT	VGDLVDLLIQ	NEFFAPASLL	LPD AVPKTAN	TLPSKEAITV	120
121	QQKQMPFCDK	DRTLMTPVQN	LEQSYMPDSD	SSPENKSLEV	SDTRFHSFSF	YELKNVTNNF	180
181	DERPISVGGN	KMGEFGFV	YKGYVNNTV	AVKKLAAMVD	ITTEELKQOF	DQEIKVMAKC	240
241	QHENVLELLG	FSSDGDDLCL	VYVYMPNGSL	LDRLSCLDGT	PPLSWHMRCK	IAQGAANGIN	300
301	FLHENHHIHR	DIKSANILLD	EAF TAKISDF	GLARASEKFA	QTVMTSRIVG	TTAYMAPEAL	360
361	RGEITPKSDI	YSFGVLEI	ITGLPAVDEH	REPQLLLDIK	EEIEDEEKT	EDYIDKMMND	420
421	ADSTSVEAMY	SVASQCLHEK	KNKRPDIKVV	QQLLQEMTAS			480

blue: IRAK4 sequence expressed in recombinant protein

¹[NCBI/Protein](https://www.ncbi.nlm.nih.gov/Protein/NP_057207.2) accession number NP_057207.2

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.