

RIPK2

receptor-interacting serine-threonine kinase 2

Recombinant Human Active Protein Kinase

HGNC Symbol: RIPK2

Synonyms: CARD3; CARDIAK; CCK; GIG30; RICK; RIP-2; RIP2

Product No.: 0676-0000-1

Lot: 003

Description: Human RIPK2, N-terminal fragment, amino acids M₁-E₂₉₉ (as in NCBI/Protein entry NP_003812.1), N-terminal GST-HIS₆ fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

Product identity: RIPK2, Lot 003, was confirmed as RIPK2 by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 63,815 Da

Expression: Baculovirus infected Sf9 cells/E.coli

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: 0.206 µg/µl

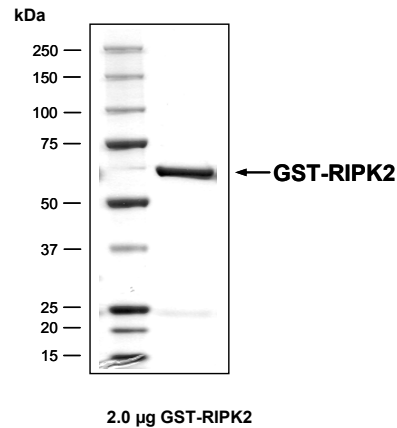
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

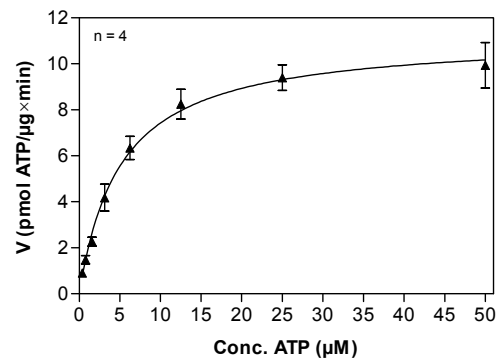
Specific kinase activity (P_i transfer): 11 pmol/µg×min

ATP-K_M: 5.1 µM

**RIPK2 Lot 003:
Coomassie stain**



**RIPK2 Lot 003:
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: RBER-CHKtide 50 µg/ml
 - Kinase: 1.0 µg/ml
- Filter binding assay
 - MSFC membrane (Millipore)

RIPK2

Product No.: 0676-0000-1

RIPK2 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQ SMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP EML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI PQID	KYLKSSKYIA	WPLQG WQATF	GGGDHPPKSD	PMGHHHHH HG	RRRASVAAGI	240
241	LVPRGSPGLD	GIYARGMNGE	AICSALPTIP	YHKLADLRYL	SRGASGTVSS	ARHADWRVQV	300
301	AVKHLHIHTP	LLDSERKDV L	REAEILHKAR	FSYILPILGI	CNEPEFLGIV	TEYMPNGSLN	360
361	ELLHRKTEYP	DVAWPLRFRI	LHEIALGVNY	LHNMTPELLH	HDLKTQNILL	DNEFHVKIAD	420
421	FGLSKWRMMS	LSQSRSSKSA	PEGGTIIYMP	PENYEPGQKS	RASIKHDIYS	YAVITWEVLS	480
481	RKQPFEDVTN	PLQIMYSVSQ	GHRPVINEES	LPYDIPHRAR	MISLIESGWA	QNPDERPSFL	540
541	KCLIELEPVL	RTFEE					600

1-218: GST **Red:** HIS6-tag **Pink:** Thrombin cleavage site **blue:** RIPK2 fragment

RIPK2 wt ¹ Amino Acid Sequence							
1	MNGEAICSAL	PTIPYHKLAD	LRYL SRGASG	TVSSARHADW	RVQVAVKHLH	IHTPLLD SER	60
61	KDVLREAEIL	HKARFSYILP	ILGICNEPEF	LGIVTEYMPN	GSLNELLHRK	TEYPDVAWPL	120
121	RFRILHEIAL	GVNYLHNMT P	PLLHDLKTQ	NILLDNEFHV	KIADFGLSKW	RMMSLSQSRS	180
181	SKSAPEGGTI	IYMPENYEP	GQKSRASIKH	DIYSYAVITW	EVLSRKQPFE	DVTNPLQIMY	240
241	SVSQGHRPVI	NEESLPYDIP	HRARMISLIE	SGWAQN PDER	PSFLKCLIEL	EPVLR T FEEI	300
301	TFLEAVIQLK	KTKLQSVSSA	IHLCDKKKME	LSLNIPVNHG	PQEESCGSSQ	LHENSGSPET	360
361	SRSLPAPQDN	DFLSRKAQDC	YFMKLHHC PG	NHSWDSTISG	SQRAAFCDHK	TTPCSSAIIN	420
421	PLSTAGNSER	LQPGIAQQWI	QSKREDIVNQ	MTEACL NQSL	DALLSRDLIM	KEDYELVSTK	480
481	PTRTSKVRQL	LDTTDIQGEE	FAKIVIVQKLK	DNKQMGLOPY	PEILVVSRS P	SLNLLQNKSM	540

blue: RIPK2 sequence expressed in fusion protein

¹NCBI/Protein accession number NP_003812.1