

RIPK4

Receptor-interacting serine-threonine kinase 4

Recombinant Human Active Protein Kinase

HGNC Symbol: RIPK4

Synonyms: ANKK2, ANKRD3, DIK, PKK, RIP4

Product No.: 1518-0000-1

Lot: 007

Description: Human RIPK4, N-terminal fragment, amino acids M₁-D₃₀₀ (as in NCBI/Protein entry NP_065690.2), N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

Product identity: RIPK4 Lot 007, was confirmed as RIPK4 by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 63,283 Da

Expression: Baculovirus infected Sf9 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

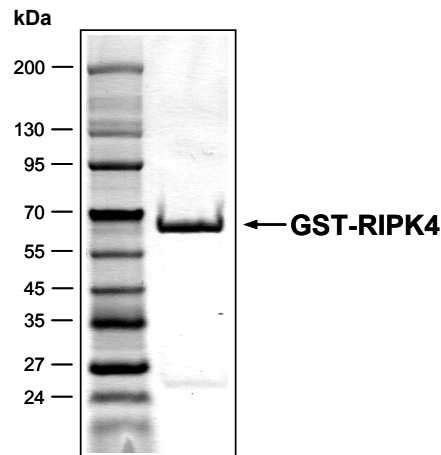
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 0.319 µg/µl (Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

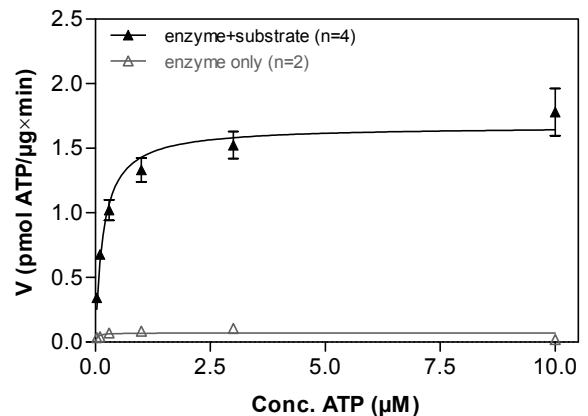
Specific kinase activity (P_i transfer): 2 pmol/µg×min
ATP-K_M: 0.17 µM

**RIPK4 Lot 007:
Coomassie stain**



2.0 µg GST-RIPK4

**RIPK4 Lot 007:
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: Casein, 20 µg / ml
 - RIPK4: 4.0 µg / ml
- Filter binding assay
 - MSFC membrane (Millipore)

Additional assay technology: RIPK4 Lot 007

was also successfully tested by ProQinase for the use with the ADP-Glo™ Kinase assay from ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details



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.

CoA V1.1 RIPK4_Lot007_V1.doc

RIPK4

Product No.: 1518-0000-1

RIPK4 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPEML	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPOID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMG HHHHHG	RDS LEVLFG	240
241	PLAMGARGRM	EGDGGTPWAL	ALLRTFDAGE	FTGWEKVGSG	GFGQVYKVRH	VHWKTWLAIK	300
301	CSPSLHVDDR	ERMELLEAK	KMEMAKFRYI	LPVYGICREP	VGLVMEYMET	GSLEKLLASE	360
361	PLPWLRFRI	IHETAVGMNF	LHCMAPPLLH	LDLKPANILL	DAHYHVKISD	FGLAKCNGLS	420
421	HSHDLSMDGL	FGTIAYLPE	RIREKSRLF	TKHDVYSFAI	VIWGVLTQKK	PFADEKNILH	480
481	IMVKVVKGHR	PELPPVCRAR	PRACSHLIRL	MQRWQGDPR	VRPTFQEITS	ETEDLCEKPD	540
541	DEVKETAHD						600

1-218: GST **Red:** HIS6-tag **Green:** 3C cleavage site **blue:** RIPK4 fragment

RIPK4 wt ¹ Amino Acid Sequence							
1	MEGDGGTPWA	LALLRTFDAG	EFTGWEKVG	GGFGQVYKVR	HVHWKTWLAI	KCSPSLHVDD	60
61	RERMELLEEA	KMEMAKFRY	ILPVYGICRE	PVGLVMEYME	TGSLEKLLAS	EPLPWLRFRI	120
121	I IHETAVGMN	FLHCMAPPLL	HLDLKPANIL	LDAHYHVKIS	DFGLAKCNGL	SHSHDLSMDG	180
181	LFGTIAYLPP	ERIREKSRLF	DTKHDVYSFA	IVIWGVLTQK	KPFADEKNIL	HIMVKVVKGH	240
241	RPELPPVCRA	RPRACSHLIR	LMQRWQGD	RVRPTFQEIT	SETEDLCEKP	DDEVKETAHD	300
301	LDVKSPPEPR	SEVVPARLKR	ASAPTFDNDY	SLSELSQLD	SGVSQAVEGP	EELSRSSSES	360
361	KLPSSGSGKR	LSGVSSVDSA	FSSRGSLSLS	FEREPSTSDL	GTTDVQKKKL	VDAIVSGDTS	420
421	KLMKILQPQD	VDLALDSGAS	LLHLAVEAGQ	EECAKWLLEN	NANPNLSNRR	GSTPLHMAVE	480
481	RRVRGVVELL	LARKISVNAK	DEDQWTALHF	AAQNGDESST	RLLEKNASV	NEVDFEGRTP	540
541	MHVACQHGE	NIVRILLRRG	VDSVSLQKDA	WLPLHYAAWQ	GHLPIVKLLA	KQPGVSVNAQ	600
601	TLDGRTPLHL	AAQRGHYRVA	RILIDLCSVD	NVCSLLAQTP	LHVAAETGHT	STARLLLRHG	660
661	AGKEAMTSDG	YTALHLAARN	GHLATVKLLV	EEKADVLARG	PLNQ TALHLA	AAHGSEVVE	720
721	ELVSADVIDL	FDEQGLSALH	LAAQRHAQT	VETLLRHGAH	INLQSLKFQG	GHGPAATLLR	780
781	RSKT						840

blue: RIPK4 sequence expressed in fusionprotein

¹NCBI/Protein accession number NP_065690.2

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.