

## ProQinase™ 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 M<sub>1</sub>-D<sub>300</sub> (as in NCBI/Protein entry NP\_065690.2), N-terminal GST-HIS<sub>6</sub> 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<sub>Fusion Protein</sub>:** 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:

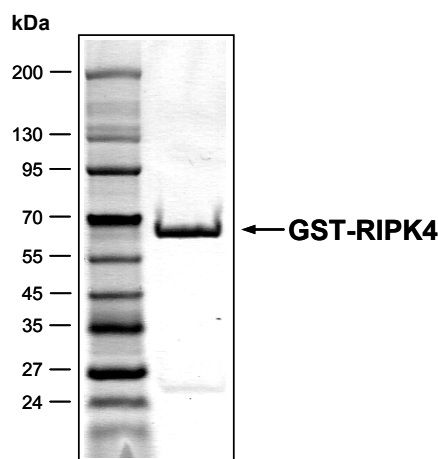
Specific kinase activity (P<sub>i</sub> transfer): 2 pmol/µg×min  
ATP-K<sub>M</sub>: 0.17 µM

**Additional assay technology:** RIPK4 Lot 007

was also successfully tested by Reaction Biology 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

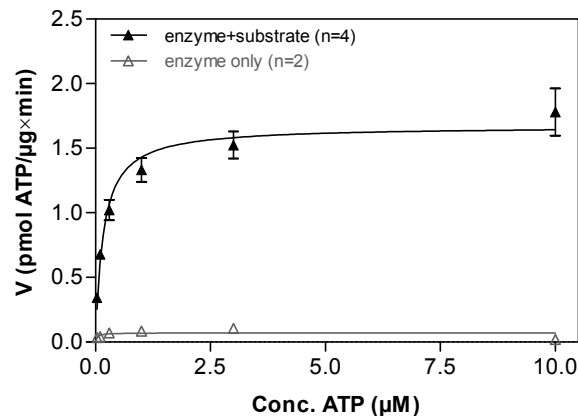


### RIPK4 Lot 007: Coomassie stain



2.0 µg GST-RIPK4

### RIPK4 Lot 007: Determination of V<sub>max</sub> and K<sub>M</sub> value for ATP



### Determination of K<sub>M</sub> value & Specific activity:

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

## ProQinase™ RIPK4

Product No.: 1518-0000-1

RIPK4 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGWYKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEML	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPOID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHG	RDSLEVLFOG	240
241	PLAMGARGRM	EGDGGTPWAL	ALLRTFDAGE	FTGWKVGSG	GFGQVYKVRH	VHWKTWLAIK	300
301	CSPSLHVDDR	ERMELLEAK	KMEMAKFRYI	LPVYGICREP	VGLVMEYMET	GSLEKLLASE	360
361	PLPWLDRFRI	IHETAVGMNF	LHCMAPPLLH	LDLKPANILL	DAHVYVKISD	FGLAKCNGLS	420
421	HSHDLSMDGL	FGTIAYLPPE	RIREKSRLF	TKHDVYSFAI	VIWGVLTQKK	PFADKCNILH	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 <sup>1</sup> Amino Acid Sequence							
1	MEGDGGTPWA	LALLRTFDAG	EFTGWKVGSG	GGFGQVYKVR	HVHWKTWLAI	KCSPSLHVDD	60
61	RERMELLEEA	KMEMAKFRY	ILPVYGICRE	PVGLVMEYME	TGSLEKLLAS	EPLPWLDRFR	120
121	IIHETAVGMN	FLHCMAPPLL	HLDLKPANIL	LDAHVYVKIS	DFGLAKCNGL	SHSHDLSMDG	180
181	LFGTIAYLPP	ERIREKSRLF	DTKHDVYSFA	IVIWGVLTQK	KPFADKCNIL	HIMVKVVKGH	240
241	RPELPPVCRA	RPRACSHLIR	LMQRWQGDPR	RVRPTFQEIT	SETEDLCEKP	DDEVKETAHD	300
301	LDVKSPPEPR	SEVVPARLKR	ASAPTFDNDY	SLSELLSQLD	SGVSQAVEGP	EELSRSSSES	360
361	KLPSSGSGKR	LSGVSSVDSA	FSSRGSLSL	FEREPSTSDL	GTTDVQKKKL	VDAIVSGDTS	420
421	KLMKILQPQD	VDLALDSGAS	LLHLAVEAGQ	EECAKWLLEN	NANPNLSNRR	GSTPLHMAVE	480
481	RRVRGVVELL	LARKISVNAK	DEDQWTALHF	AAQNGDESST	RLLLEKNASV	NEVDFEGRTP	540
541	MHVACQHQQE	NIVRIILLRRG	VDVSLQKDA	WLPLHYAAWQ	GHLPIVKLLA	KQPGVSVNAQ	600
601	TLDGRTPLHL	AAQRGHYRVA	RILIDLCSVD	NVCSLLAQTP	LHVAAETGHT	STARLLLHRG	660
661	AGKEAMTSDG	YTALHLAARN	GHLATVKLLV	EEKADVLARG	PLNQ TALHLA	AAHGHSEVVE	720
721	ELVSADVIDL	FDEQGLSALH	LAAQGRHAQT	VETLLRHGAH	INLQSLKFQG	GHGPAATLLR	780
781	RSKT						840

**blue:** RIPK4 sequence expressed in fusionprotein

<sup>1</sup>NCBI/Protein accession number NP\_065690.2

**Recombinant Proteins**