

## ProQinase™ PAK4

p21 protein (Cdc42/Rac)-activated kinase 4

Recombinant Human Active Protein Kinase

HGNC Symbol: PAK4

Synonyms: n/a

Product No.: 0365-0000-1

Lot: 004

**Description:** Human PAK4, full length, amino acids M<sub>1</sub>-R<sub>591</sub> (as in NCBI/Protein entry NP\_005875.1), N-terminal GST-HIS<sub>6</sub> fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

**Product identity:** PAK4 Lot 004, was confirmed as PAK4 by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW<sub>Fusion Protein</sub>:** 93,968 Da

**Expression:** Baculovirus infected Sf9 cells

**Purification:** GST-Affinity Chromatography

**Activation:** This kinase was not activated by special procedures

**Storage buffer:** 50 mM Tris-HCl pH 8.0, 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:** 0.154 µg/µl

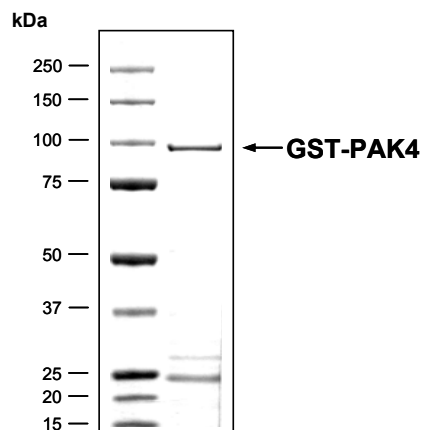
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

**Biochemical Parameters:**

Specific kinase activity (P<sub>i</sub> transfer): 3 pmol/µg×min

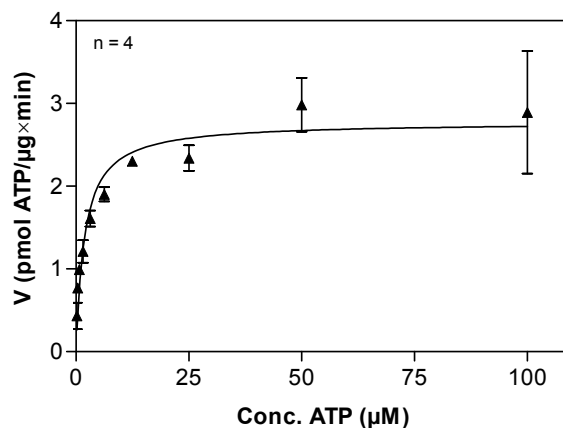
ATP-K<sub>M</sub>: 1.9 µM

### PAK4 Lot 004: Coomassie stain



2.0 µg GST-PAK4

### PAK4 Lot 004: 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: tetra(LRRWSLG), 50 µg/ml
  - PAK4: 4.0 µg/ml
- Filter binding assay
  - MSPH membrane (Millipore)

## ProQinase™ PAK4

Product No.: 0365-0000-1

PAK4 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPEML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPOID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMCHHHHHHG	RRRASVAAGI	240
241	LVPRGSPGLD	GIYARGIQAS	MFGKRKKRVE	ISAPSNFEHR	VHTGFDQHEQ	KFTGLPRQWQ	300
301	SLIEESARRP	KPLVDPACIT	SIQPGAPKTI	VRGSKGAKDG	ALTLILLDEFE	NMSVTRSNSL	360
361	RRDSPPPPAR	ARQENGMPEE	PATTARGGPG	KAGSRGRFAG	HSEAGGGSGD	RRRAGPEKRP	420
421	KSSREGSGGP	QESSRDKRPL	SGPDVGTQPQ	AGLASGAKLA	AGRPFNTYPR	ADTDHPSRGA	480
481	QGEPHDVAPN	GPSAGGLAIP	QSSSSSRPP	TRARGAPSPG	VLGPHASEPQ	LAPPACTPAA	540
541	PAVPGPPGPR	SPQREPQRVS	HEQFRAALQL	VDPGDPRS	LDNFIKIGEG	STGIVCIATV	600
601	RSSGKLVAVK	KMDLRKQRR	ELLFNEVVIM	RDYQHENVVE	MYSYLVGDE	LWVMEFLEG	660
661	GALTDIVTHT	RMNEEQIAAV	CLAVLQALSV	LHAQGVHRD	IKSDSILLTH	DGRVKLSDFG	720
721	FCAQVSKEVP	RRKSLVGTPY	WMAPELISRL	PYGPEVDIWS	LGIMVIEMVD	GEPPYFNEPP	780
781	LKAMKIRDN	LPPRLKNLHK	VSPSLKGFLD	RLLRDPAQR	ATAAELLKHP	FLAKAGPPAS	840
841	IVPLMRQNR	R					900

1-218: GST Red: HIS6-tag Pink: Thrombin cleavage site blue: PAK4

PAK4 wt <sup>1</sup> Amino Acid Sequence							
1	MFGKRKKRVE	ISAPSNFEHR	VHTGFDQHEQ	KFTGLPRQWQ	SLIEESARRP	KPLVDPACIT	60
61	SIQPGAPKTI	VRGSKGAKDG	ALTLILLDEFE	NMSVTRSNSL	RRDSPPPPAR	ARQENGMPEE	120
121	PATTARGGPG	KAGSRGRFAG	HSEAGGGSGD	RRRAGPEKRP	KSSREGSGGP	QESSRDKRPL	180
181	SGPDVGTQPQ	AGLASGAKLA	AGRPFNTYPR	ADTDHPSRGA	QGEPHDVAPN	GPSAGGLAIP	240
241	QSSSSSRPP	TRARGAPSPG	VLGPHASEPQ	LAPPACTPAA	PAVPGPPGPR	SPQREPQRVS	300
301	HEQFRAALQL	VDPGDPRS	LDNFIKIGEG	STGIVCIATV	RSSGKLVAVK	KMDLRKQRR	360
361	ELLFNEVVIM	RDYQHENVVE	MYSYLVGDE	LWVMEFLEG	GALTDIVTHT	RMNEEQIAAV	420
421	CLAVLQALSV	LHAQGVHRD	IKSDSILLTH	DGRVKLSDFG	FCAQVSKEVP	RRKSLVGTPY	480
481	WMAPELISRL	PYGPEVDIWS	LGIMVIEMVD	GEPPYFNEPP	LKAMKIRDN	LPPRLKNLHK	540
541	VSPSLKGFLD	RLLRDPAQR	ATAAELLKHP	FLAKAGPPAS	IVPLMRQNR	R	600

blue: PAK4 sequence expressed in fusionprotein

<sup>1</sup>NCBI/Protein accession number NP\_005875.1

Recombinant Proteins