

ProQinaseTM GAK

cyclin G associated kinase

Recombinant Human Active Protein Kinase

HGNC Symbol: GAK

Synonyms: Auxilin-2, DNAJC26

Product No.: 1974-0000-1

Lot: 003

Description: Human Kinase, N-terminal fragment, amino acids M₁-V₄₀₀ (as in [NCBI/Protein](#) entry NP_005246.2), untagged, expressed in Sf9 insect cells

Product identity: GAK Lot 003, has been verified by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 44510 Da

Expression host: Sf9 insect cells

Purification: GST-Affinity Chromatography

Activation: in vitro auto activation

Storage buffer: 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 20 % glycerol

Storage temperature: -80°C

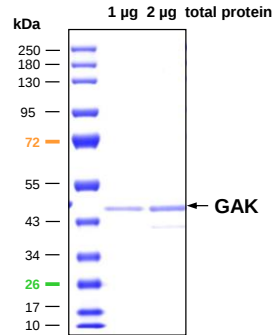
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10 µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 0.466 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

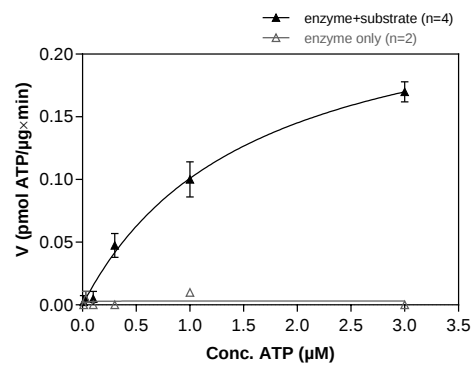
Biochemical Parameters:

Specific kinase activity (P_i transfer): 0.3 pmol/µg*min
ATP-K_M: 1.6 µM

**GAK Lot 003:
Coomassie stain**



**GAK Lot 003:
Determination of V_{max} and K_M value for ATP**



- 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: ERK2-K54R 80 µg/ml
 - Kinase: 4 µg/ml

Assay technology:
Radiometric filter binding assay
MSFC membrane (96 well plate, Millipore)

Sequence information

GAK Recombinant Fusion Protein Amino Acid Sequence

1	GPMSLLQSAL	DFLAGPGSLG	GASGRDQSDF	VGQTVELGEL	RLRVRRVLAE	GGFAFVYEAQ	60
61	DVGSGREYAL	KRLLSNEEEK	NRAIIQEVCF	MKKLSGHPNI	VQFCSAASIG	KEESDTGQAE	120
121	FLLLTELCKG	QLVEFLKKME	SRGPLSCDTV	LKIFYQTCRA	VQHMHRQKPP	IIHRDLKVEN	180
181	LLLSNQGTIK	LCDFGSATTI	SHYPDYSWSA	QRRALVEEEI	TRNTTPMYRT	PEIIDLYSNF	240
241	PIGEKQDIWA	LGCILYLLCF	RQHPFEDGAK	LRIVNGKYSI	PPHDTQYTVF	HSLIRAMLQV	300
301	NPEERLSIAE	VVHQLQEIAA	ARNVNPKSPI	TELEQNGGY	GSATLSRGPP	PPVGAGSGY	360
361	SGGLALAEYD	PYGGFLDIL	RGGTERLFTN	LKDTSSKVIQ	SV		420

1-2: leagcy of tag-cleavage blue: GAK fragment

GAK wt¹ Amino Acid Sequence

1	MSLLQSALDF	LAGPGSLGGA	SGRDQSDFVG	QTVELGELRL	RVRRLAEGG	FAFVYEAQDV	60
61	SGGREYALKR	LLSNEEEKNR	AIIEVCFMK	KLSGHPNIVQ	FCSAASIGKE	ESDTGQAEFL	120
121	LLTELCKGQL	VEFLKKMESR	GPLSCDTVLK	IFYQTCRAVQ	HMHROKPPII	HRDLKVENLL	180
181	LSNQGTIKLC	DFGSATTISH	YPDYSWSAQR	RALVEEEITR	NTPPMYRTPE	IIDLYSNFPI	240
241	GEKQDIWALG	CILYLLCFRQ	HPFEDGAKLR	IVNGKYSIPP	HDTQYTVFHS	LIRAMLQVNP	300
301	EERLSIAEVV	HQQLFIAAAR	NVNPKSPITE	LLEQNGGYGS	ATLSRGPPPP	VGPAGSGYSG	360
361	GLALAEYDQP	YGGFELDILRG	GTERLFTNLK	DTSSKVIQSV	ANYAKGDLDI	SYITSRIAVM	420
421	SFPAEGVESA	LKNNIEDVRL	FLDSKHPGHY	AVYNLSPRTY	RPSRFHNRVS	ECGWAARRAP	480
481	HLHTLYNICR	NMHAWLRQDH	KNVCVVHCMD	GRAASAVAVC	SFLCFCRLFS	TAAEAAYMFS	540
541	MKRCPPGIWP	SHKRYIEYMC	DMVAEEPITP	HSKPILVRVAV	VMTVPVPLFSK	QRSGCRPFCE	600
601	VYVGDERVAS	TSQEYDKMRD	FKIEDGKAVI	PLGVTVQGDV	LIVIIYHARST	LGGRLLQAKMA	660
661	SMKMFQIQFH	TGFVPRNATT	VKFAKYDLDA	CDIQEKYPDL	FQVNLEVEVE	PRDRPSREAP	720
721	PWENSSMRGL	NPKILFSSRE	EQQDILSKFG	KPELRPQSGS	TAQYDAGAGS	PEAEPTDSDS	780
781	PPSSADASRL	FLHTLDWQEE	KEATEGAENA	SSKESSESALM	EDRDESEVSD	EGGSPISSEG	840
841	QEPRADPEPP	GLAAGLVQQD	LVFEVETPAV	LPEVPVQEDG	VDLLGLHSEV	GAGPAVPPQA	900
901	CKAPSSNTDL	LSCLLGPEEA	ASQGPPEDLL	SEDPDLLLASP	APPLSVQSTP	RGGPPAAADP	960
961	FGPLLPSSGN	NSQPCSNPDL	FGEFLNSDSV	TVPPSFPSAH	SAPPPSCSAD	FLHLGDLPGE	1020
1021	PSKMTASSSN	PDLLGGWAAW	TETAASAVAP	TPATEGPLFS	PGGPAPCGS	QASWTKSQNP	1080
1081	DPFADLGDL	SLQGGSPAGF	PPGGFIPKTA	TTPKGSSSWQ	TSRPPAQGAS	WPPQAKPPPK	1140
1141	ACTQPRPNYA	SNFSVIGARE	ERGVRAPSGA	QPKVSENDF	EDLLSNQGSF	SRSDKKGPKT	1200
1201	IAEMRKQDLA	KDTPDLKLKL	LDWIEGKERN	IRALLSTLHT	VLWDGESRWT	PVGMADLVAP	1260
1261	EQVKKHYRRA	VLAHPDKAA	GQPYEQHAKM	IFMELNDAWS	EFENOGSRPL	F	1320

blue: GAK sequence expressed in recombinant protein

¹NCBI/Protein accession number NP_005246.2