

ProQinase™ KIT V560G

KIT proto-oncogene, receptor tyrosine kinase

Recombinant Human Active Protein Kinase

HGNC Symbol: KIT

Synonyms: CD117, PBT, SCFR, c-Kit

Product No.: 1046-0000-1

Lot: 002

Description: Human KIT, C-terminal fragment, amino acids T₅₄₄-V₉₇₆ (as in [NCBI/Protein](#) entry NP_000213.1) with a V560G mutation, N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

Product identity: KIT V560G Lot 002 was confirmed as KIT by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 77,415 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, 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: 0.101 µg/µl

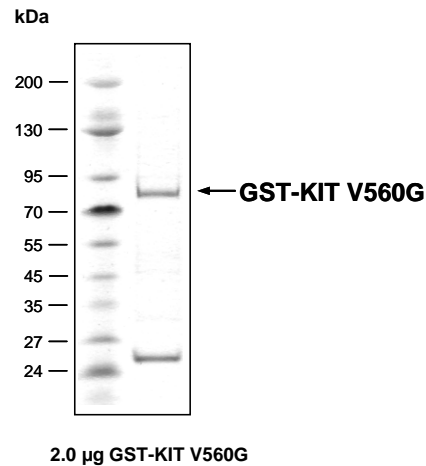
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

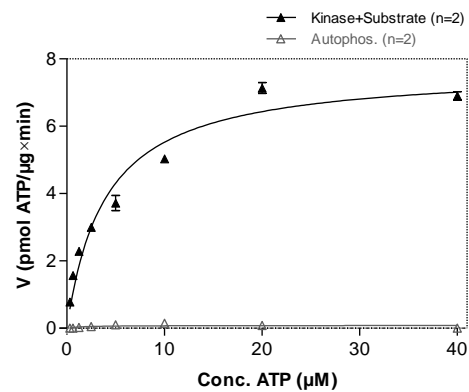
Specific kinase activity (P_i transfer): 8 pmol/µg × min

ATP-K_M: 3.9 µM

KIT V560G Lot 002: Coomassie stain



KIT V560G Lot 002: 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: TRK-C derived peptide 80 µg/ml
Kinase: 4 µg/ml
- Filter binding assay
MSPH membrane (Millipore)

Additional assay technology:

KIT V560G Lot 002 was also successfully tested by Reaction Biology for the use with the ADP-Glo™ Kinase assay from Promega ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details

ProQinase™ KIT V560G

Product No.: 1046-0000-1

GST-KIT V560G Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPIQID	KYLKSSKYIA	WPLQGWAQTF	GGGDHPPKSD	PMG HHHHHG	RDS LEVLFQG	240
241	PLAMGTYKYL	QKPMYEVQWK	V E E I N G N N Y	V I D P T Q L P Y	D H K W E F P R N R	L S F G K T L G A G	300
301	A F G K V E A T	Y G L I K S D A A M	T V A V K M L K P S	A H L T E R E A L M	S E L K V L S Y L G	N H M N I V N L L G	360
361	A C T I G G P T L V	I T E Y C C Y G D L	L N F L R R K R D S	F I C S K Q E D H A	E A A L Y K N L L H	S K E S S C S D S T	420
421	N E Y M D M K P G V	S Y V V P T K A D K	R R S V R I G S Y I	E R D V T P A I M E	D D E L A L D L E D	L L S F S Y Q V A K	480
481	G M A F L A S K N C	I H R D L A A R N I	L L T H G R I T K I	C D F L A R D I K	N D S N Y V V K G N	A R L P V K W M A P	540
541	E S I F N C V Y T F	E S D V W S Y G I F	L W E L F S L G S S	P Y P G M P V D S K	F Y K M I K E G F R	M L S P E H A P A E	600
601	M Y D I M K T C W D	A D P L K R P T F K	Q I V Q L I E K Q I	S E S T N H I Y S N	L A N C S P N R Q K	P V V D H S V R I N	660
661	S V G S T A S S Q	P L L V H D D V					720

1-218: GST **Red**: HIS6-tag **Green**: 3C cleavage site **blue**: KIT fragment **boxed**: V560G

KIT wt ¹ Amino Acid Sequence							
1	MARGAWDF	LCVLLLLLRV	QTGSSQPSVS	PGEPSPPSIH	PGKSDLIVRV	GDEIRLLCTD	60
61	PGFVKWTFEI	LDETENENQN	EWITEKAEAT	NTGKYTCTNK	HGLSNSIYVF	VRDPAKFLV	120
121	DRSLYGKEDN	DTLVRCPDLD	PEVTNYSKLG	CQGKPLPKDL	RFIPDPKAGI	MIKSVKRAYH	180
181	RLCLHCSVDQ	EGKSVLSEKF	ILKVRPAFKA	VPVSVSKAS	YLLREGEEFT	VTCTIKDVSS	240
241	SVYSTWKREN	SQTKLQEKYN	SWHHGDFNYE	RQATLTISSA	RVNDSGVFMC	YANNTFGSAN	300
301	VTTTLEVVDK	GFINIFPMIN	TTVFVNDGEN	VDLIVEYEAF	PKPEHQWIY	MNRTFTDKWE	360
361	DYPKSENESEN	IRYVSELHLT	RLKGTEGGTY	TFLVSNSDVN	AAIAFNIVVN	TKPEILTYDR	420
421	LVNGMLQCVA	AGFPEPTIDW	YFCPGTEQRC	SASVLPVDVQ	TLNSSGPPFG	KLVVQSSIDS	480
481	SAFKHNGTVE	CKAYNDVGKT	SAYFNFAFKG	NNKEQIHPHT	LFTPLLIGFV	IVAGMMCIIV	540
541	MILTYKYLQK	P M Y E V Q W K V	E E I N G N N Y V Y	I D P T Q L P Y D H	K W E F P R N R L S	F G K T L G A G A F	600
600	G K V E A T A Y G	L I K S D A A M T V	A V K M L K P S A H	L T E R E A L M S E	L K V L S Y L G N H	M N I V N L L G A C	660
661	T I G G P T L V I T	E Y C C Y G D L L N	F L R R K R D S F I	C S K Q E D H A E A	A L Y K N L L H S K	E S S C S D S T N E	720
721	Y M D M K P G V S Y	V V P T K A D K R R	S V R I G S Y I E R	D V T P A I M E D D	E L A L D L E D L L	S F S Y Q V A K G M	780
781	A F L A S K N C I H	R D L A A R N I L	T H G R I T K I C D	F L G A R D I K N D	S N Y V V K G N A R	L P V K W M A P E S	840
841	I F N C V Y T F E S	D V W S Y G I F L W	E L F S L G S S P Y	P G M P V D S K F Y	K M I K E G F R M L	S P E H A P A E M Y	900
901	D I M K T C W D A D	P L K R P T F K Q I	V Q L I E K Q I S E	S T N H I Y S N L A	N C S P N R Q K P V	V D H S V R I N S V	960
961	G S T A S S Q P L	L V H D D V					1020

blue: KIT sequence expressed in recombinant protein **Red**: variant in recombinant protein

¹[NCBI/Protein](#) accession number NP_000213.1