

## ProQinase™ MET L1195V

met proto-oncogene

Recombinant Human Active Protein Kinase

HGNC Symbol: MET

Synonyms: c-MET, HGFR

Product No.: 1651-0000-1

Lot: 001

**Description:** Human MET C-terminal fragment, amino acids K<sub>956</sub>-S<sub>1390</sub> (as in [NCBI/Protein](#) entry NP\_000236.2), L<sub>1195</sub>V point mutant, N-terminal GST-HIS<sub>6</sub> fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

**Product identity:** MET L1195V Lot 001 product identity was confirmed by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW<sub>Fusion Protein</sub>:** 78,773 Da

**Expression host:** Sf9 insect 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:** 0.173 µg/µl

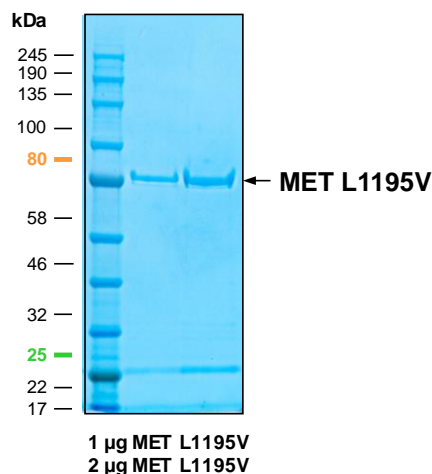
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

**Biochemical Parameters:**

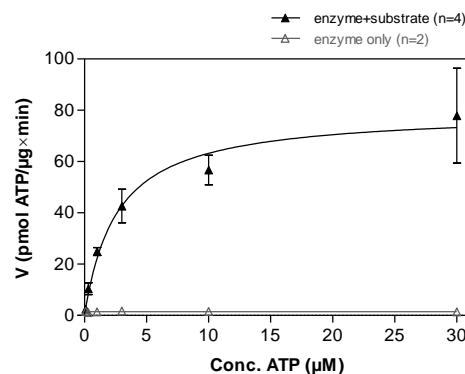
Specific kinase activity (P<sub>i</sub> transfer): 80 pmol/µg × min

ATP-K<sub>M</sub>: 2.6 µM

**MET L1195V Lot 001:**  
Coomassie stain



**MET L1195V Lot 001:**  
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: TRK-C derived peptide 20 µg/ml

Kinase: 1 µg/ml

• Filter binding assay

MSPH membrane (Millipore)

**Additional assay technology:**

MET L1195V Lot 001 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™ MET L1195V

Product No.: 1651-0000-1

GST-MET L1195V Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQDMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMG <b>HHHHHG</b>	RRRASVAAGI	240
241	<b>LVPRG</b> SPGLD	GICSIIEFK	<b>RKQIKDLGSE</b>	<b>LVRVDARVHT</b>	<b>PHLDRLVSAR</b>	<b>SVSPTTEMVS</b>	300
301	<b>NESVDYRATF</b>	<b>PEDQFPNSSQ</b>	<b>NGSCRQVQYP</b>	<b>LTDMSPI LTS</b>	<b>GSDISSPLL</b>	<b>QNTVHIDL SA</b>	360
361	<b>LNPELVQAVQ</b>	<b>HVVIGPSSLI</b>	<b>VHFNEVIGRG</b>	<b>HFGCVYHGTL</b>	<b>LDNDGKKIHC</b>	<b>AVKSLNRTD</b>	420
421	<b>IGEVSQFLTE</b>	<b>GIIMKDFSHP</b>	<b>NVLSLLGICL</b>	<b>RSEGSPLVVL</b>	<b>PYMKHGDLRN</b>	<b>FIRNETHNPT</b>	480
481	<b>VKDLIGFGLQ</b>	<b>VAKGMKYVAS</b>	<b>KKFVHRDLAA</b>	<b>RNCMLDEKFT</b>	<b>VKVADFGLAR</b>	<b>DMYDKEYYSV</b>	540
541	<b>HNKTKAKLPV</b>	<b>KWMALESLOT</b>	<b>QKFTTKSDVW</b>	<b>SFGVLLWELM</b>	<b>TRGAPPYDPV</b>	<b>NTFDITVYLL</b>	600
601	<b>QGRRLLOPEY</b>	<b>CPDPLYEVML</b>	<b>KCWHPKAEMR</b>	<b>PSFSELVSRI</b>	<b>SAIFSTFIGE</b>	<b>HYVHV NATYV</b>	660
661	<b>NVKCVAPYPS</b>	<b>LLSSEDNADD</b>	<b>EVDTRPASFW</b>	<b>ETS</b>			720

1-218: GST **Red**: HIS6-tag **Pink**: Thrombin cleavage site **blue**: MET fragment **boxed**: L1195V point mutation

MET wt <sup>1</sup> Amino Acid Sequence							
1	MKAPAVLAPG	ILVLLFTLVQ	RSNGECKEAL	AKSEMNVNMK	YQLPNFTAET	PIQNVILHEH	60
61	HIFLGATNYI	YVLNEEDLQK	VAEYKTGPVL	EHPDCFPQD	CSSKANLSGG	VWKDINIMMAL	120
121	VVDYYDDQL	ISCGSVNRGT	CQRHVFPNH	TADIQSEVHC	IFSPQIEEPS	QCPDCVVSAL	180
181	GAKVLSVKD	RFINFFVGN	INSSYFPDHP	LHSISVRLK	ETKDGFMFLT	DQSYIDLPE	240
241	FRDSYPIKYV	HAFESNNFY	FLTVQRETL	AQTFHTRIR	FCSINSLHS	YMEMPLECIL	300
301	TEKRKKRSTK	KEVFNILQAA	YVSKPGAQLA	RQIGASLND	ILFGVFAQSK	PDSAEPMDRS	360
361	AMCAFPKIYV	NDFFNKIVNK	NNVRCLQHFY	GNHEHCENR	TLLRNSSGCE	ARRDEYRTEF	420
421	TTALQRVDF	MGQFSEVLLT	SISTFIKGD	TIANLGTSEG	RFMQVVVSR	GPSTPHVNF	480
481	LDSHPVSPEV	IVEHTLNQNG	YTLVITGKI	TKIPLNGLC	RHFQSCSQCL	SAPPFVQCGW	540
541	CHDKCVRSEE	CLSGTWTQOI	CLPAIYKVP	NSAPLEGGTR	LTICGWDFGF	RRNNKFDLKK	600
600	TRVLLGNESC	TLTLESTMN	TLKCTVGPAM	NKHFNMSII	SNGHGTQYS	TFSYVDPVIT	660
661	SISPKYGPMA	GGTLLTLTGN	YLNNGNSRHI	SIGGKTCTLK	SVSNSILECY	TPAQTISTEF	720
721	AVKCLKIDLAN	RETSIFSIRE	DPIVYIEHPT	KSFISGGSTI	TGVGKNLSV	SVPRMIVNH	780
781	EAGRNFVAC	QHRNSSEIIC	CTTPSLQQLN	LQLPLKTKAF	FMLDGILSKY	FDLIYVHNPV	840
841	FKPFKPVMI	SMGNENVLEI	KGNDIDPEAV	KGEVLKVGNK	SCENIHLHSE	AVLCTVPNL	900
901	LKLNSELNIE	WKQAISSTVL	GKVIVQPDQN	FTGLIAGVVS	ISTALLLLLG	FFLWL <b>KRRQ</b>	960
961	<b>IKDLGSELVR</b>	<b>YDARVHTPHL</b>	<b>DRLVSARSVS</b>	<b>PTEMVSNES</b>	<b>VDYRATFPED</b>	<b>QFPNSSQNGS</b>	1020
1021	<b>CRQVQYPLTD</b>	<b>MSPILTSGDS</b>	<b>DISSPLLQNT</b>	<b>VHIDL SALNP</b>	<b>ELVQAVQHVV</b>	<b>IGPSSLI VHF</b>	1080
1081	<b>NEVIGRGHFG</b>	<b>CVYHGTL LDN</b>	<b>DGKKIHC AVK</b>	<b>SLNRITDIGE</b>	<b>VSQFLTEGII</b>	<b>MKDFSHPNVL</b>	1140
1141	<b>SLLGICLRSE</b>	<b>GSPLVLPYM</b>	<b>KHGDLRNFIR</b>	<b>NETHNPTVKD</b>	<b>LIGFGLQVAK</b>	<b>GMYLASKKF</b>	1200
1201	<b>VHRDLAARNC</b>	<b>MLDEKFTVKV</b>	<b>ADFGLARDMY</b>	<b>DKEYYSVHNK</b>	<b>TGAKLPVKWM</b>	<b>ALESLOTQKF</b>	1260
1261	<b>TTKSDVWSFG</b>	<b>VLLWELMTRG</b>	<b>APPYDPVNTF</b>	<b>DITVYLLQGR</b>	<b>RLLOPEYCPD</b>	<b>PLYEVMLKCW</b>	1320
1321	<b>HPKAEMRPSF</b>	<b>SELVSRISAI</b>	<b>FSTFIGEHYV</b>	<b>HVNATYVNVK</b>	<b>CVAPYPSLLS</b>	<b>SEDNADDEVD</b>	1380
1381	<b>TRPASFWETS</b>						1440

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

<sup>1</sup>[NCBI/Protein](#) accession number NP\_000236.2

**Please notice:**

Variant amino acid numbering beginning with Ser755 when referring to GenBank accession number J02958 / [NCBI/Protein](#) accession number AAA59591.1 (additional 18 aa exon between S755/G756, frequently found in the literature)