

## ProQinase™ MET M1250T

met proto-oncogene

Recombinant Human Active Protein Kinase

HGNC Symbol: MET

Synonyms: c-MET, HGFR

Product No.: 0982-0000-1

Lot: 002

**Description:** Human MET C-terminal fragment, amino acids K<sub>956</sub>-S<sub>1390</sub> (as in NCBI/Protein entry NP\_000236.2), M1250T mutant, N-terminal GST-HIS<sub>6</sub> fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

**Product identity:** MET M1250T Lot 002 product identity was confirmed by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW<sub>Fusion Protein</sub>:** 78,757 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:** 0.504 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

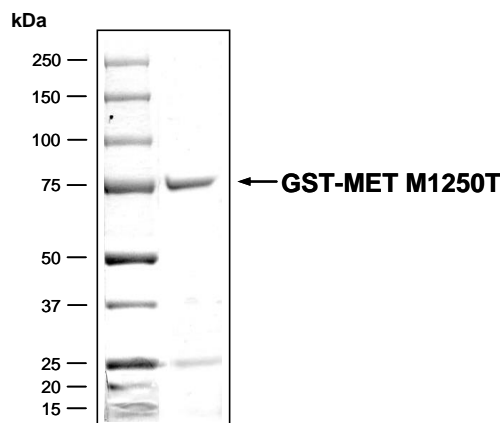
**Biochemical Parameters:**

Specific kinase activity (P<sub>i</sub> transfer): 97 pmol/µg x min  
ATP-K<sub>M</sub>: 1.4 µM

**Additional assay technology:** MET M1250T Lot 002 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

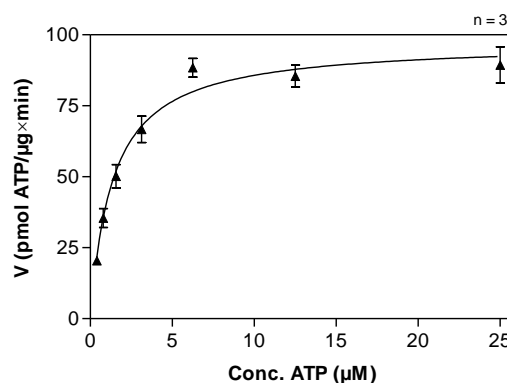


**MET M1250T Lot 002:  
Coomassie stain**



2.0 µg GST-MET M1250T

**MET M1250T Lot 002:  
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: 1.0 µg/ml
- Filter binding assay
  - MSPH membrane (Millipore)

Recombinant Proteins

# ProQinase™ MET M1250T

Product No.: 0982-0000-1

Recombinant Proteins

| MET M1250T 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 | WPLQGWQATF | GGGDHPPKSD  | PMGHHHHHG   | RRRASVAAGI | 240 |
| 241   | LVPRGSPGLD  | GICSIEEFKK | RKQIKDLGSE | LVRYDARVHT  | PHLDRLVSAR  | SVSPTEMVSV | 300 |
| 301   | NESVDYRATF  | PEDQFPNSSQ | NGSCRQVQYP | LTDMSPIILTS | GSDDISSPLL  | QNTVHIDLSA | 360 |
| 361   | LNPELVQAVQ  | HVVIGPSSLI | VHFNEVIGRG | HFGCVYHGTL  | LDNDGKKIHC  | AVKSLNRITD | 420 |
| 421   | IGEVSQFLTE  | GIIMKDFSHP | NVLSLLGICL | RSEGSPLVVL  | PYMKHGDLRN  | FIRNETHNPT | 480 |
| 481   | VKDLIGFGLQ  | VAKGMKYLAS | KKFVHRDLAA | RNCMLDEKFT  | VKVADEFGLAR | DMYDKEYYSV | 540 |
| 541   | HNKRTGAKLPV | KWTALESLOT | QKFTTKSDVM | SFGVLLWELM  | TRGAPPYPDV  | NTFDITVYLL | 600 |
| 600   | QGRRLAQPEY  | CPDPLYEVM  | KCWHPKAEMR | PSFSELVSRI  | SAIFSTFIGE  | HYVHVNATYV | 660 |
| 661   | NVKCVAPYPS  | LLSSEDNADD | EVDTRPASFW | ETS         |             |            | 720 |

1-218: GST    Red: HIS6-tag    Pink: Thrombin cleavage site    blue: MET fragment    boxed: M1250T

| MET wt <sup>1</sup> Amino Acid Sequence |             |             |            |            |            |            |      |
|---|-------------|-------------|------------|------------|------------|------------|------|
| 1                                       | MKAPAVLAPG  | ILVLLFTLVQ  | RSNGECKEAL | AKSEMNVNMK | YQLPNFTAET | PIQNVILHEH | 60   |
| 61                                      | HIFLGATNYI  | YVLNEEDLQK  | VAEYKTGPVL | EHPDCFCQD  | CSSKANLSGG | VWVDNINMAL | 120  |
| 121                                     | VVDYYDDQL   | ISCGSVNRGT  | CQRHVFPHNH | TADIQSEVHC | IFSPQIEEPS | QCPDCVVSAL | 180  |
| 181                                     | GAKVLSSVKD  | RFINFFVGN   | INSSYFPDHP | LHSISVRRLK | ETKDGFMFLT | DQSYIDVLPE | 240  |
| 241                                     | FRDSYPIKYV  | HAFESNNFIY  | FLTVQRETLD | AQTFHTRIIR | FCSINSGLHS | YMEMPLECIL | 300  |
| 301                                     | TEKRKRKSTK  | KEVFNILQAA  | YVSKPGAQLA | RQIGASLND  | ILFGVFAQSK | PDSAEPMDRS | 360  |
| 361                                     | AMCAFPKIVY  | NDFFNKIVNK  | NNVRCLQHFY | GNHEHCENR  | TLLRNSGCE  | ARRDEYRTEF | 420  |
| 421                                     | TTALQRVDLF  | MGQFSEVLLT  | SISTFIKQDL | TIANLGTSEG | RHFQSCSQCL | SAPPFVQCGW | 480  |
| 481                                     | LDSHPVSPEV  | IVEHTLNQNG  | YTLVITGKKI | TKIPLNGLGC | RHFQSCSQCL | SAPPFVQCGW | 540  |
| 541                                     | CHDKCVRSEE  | CLSGTWTQQI  | CLPAIYKVPF | NSAPLEGGTR | LTICGWDFGF | RRNNKFDLKK | 600  |
| 600                                     | TRVLLGNESC  | TLTLSESTMN  | TLKCTVGPAM | NKHFNMSIII | SNGHGTQYS  | TFSYVDPVIT | 660  |
| 661                                     | SISPKYGPMA  | GGTLLTLTGN  | YLNNGNSRHI | SIGGKTCTLK | SVSNSILECY | TPAQTISTEF | 720  |
| 721                                     | AVKCLKIDLAN | RETSIFSIRE  | DPIVYIEIHT | KSFISGGSTI | TGVGKNLNSV | SVPRMVINVH | 780  |
| 781                                     | EAGRNFVAC   | QHRNSSEIIC  | CTTPSLQQLN | LQLPLKTKAF | FMLDGILSKY | FDLIYVHNPV | 840  |
| 841                                     | FKPFKPVMI   | SMGNENVLEI  | KGNDIDPEAV | KGEVLKVGNK | SCENIHLHSE | AVLCTVPNDL | 900  |
| 901                                     | LKLNSELNIE  | WKQAISSTVL  | GKVIVQPDQN | FTGLIAGVVS | ISTALLLLL  | FFLWLKRRKQ | 960  |
| 961                                     | IKDLGSELVR  | YDARVHTPHL  | DRLVSARSVS | PTTEMVSNES | VDYRATFPED | QFPNSSQNGS | 1020 |
| 1021                                    | CRQVQPLTD   | MSPILTSQDS  | DISSPLLQNT | VHIDLSALNP | ELVQAVQHV  | IGPSSLIVHF | 1080 |
| 1081                                    | NEVIGRGHFG  | CVYHGTLN    | DGKKIHCAVK | SLNRITDIGE | VSQFLTEGII | MKDFSHPNVL | 1140 |
| 1141                                    | SLLGICLRSE  | GSPLVLPYM   | KHGDLRNFIR | NETHNPTVKD | LIGFGLQVAK | GMKYLASKKF | 1200 |
| 1201                                    | VHRDLAARNC  | MLDEKFTVKV  | ADFGLARDMY | DKEYYSVHNK | TGAKLPVKWM | ALESLOTQKF | 1260 |
| 1261                                    | TTKSDVWSFG  | VLLWELMTRG  | APPYPDVNTF | DITVYLLQGR | RLLOPEYCPD | PLYEVMLKCW | 1320 |
| 1321                                    | HPKAEMRPSF  | SELVSRIASAI | FSTFIGEYHV | HVNATYVNVK | CVAPYPSLLS | SEDNADDEVD | 1380 |
| 1381                                    | TRPASFWETS  |             |            |            |            |            | 1440 |

blue: MET sequence expressed in fusion protein    Red: variant in fusion 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 J02958 (additional 18 aa exon between S755/G756, frequently found in the literature)