

## ProKinase™ MET Y1230H

met proto-oncogene

Recombinant Human Active Protein Kinase

HGNC Symbol: MET

Synonyms: c-MET, HGFR

Product No.: 0976-0000-1

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), Y1230H mutant, N-terminal GST-HIS<sub>6</sub> fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

**Product identity:** MET Y1230H Lot 001 product identity was confirmed by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW**<sub>Fusion Protein</sub>: 78,761 Da

**Expression:** Baculovirus infected Sf9 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:** 0.295 µg/µl

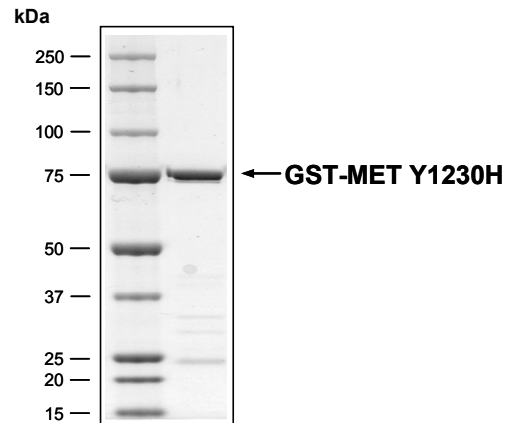
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

**Biochemical Parameters:**

Specific kinase activity (P<sub>i</sub> transfer): 116 pmol/µg × min

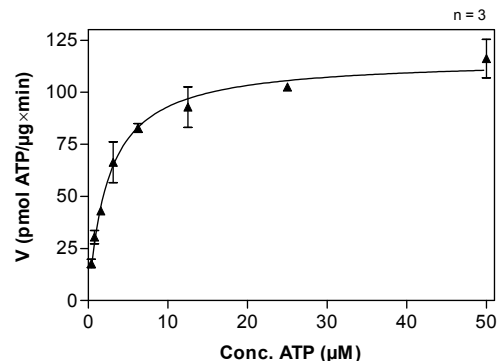
ATP-K<sub>M</sub>: 2.5 µM

**MET Y1230H Lot 001:  
Coomassie stain**



2.0 µg GST-MET Y1230H

**MET Y1230H 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: 1.0 µg/ml
- Filter binding assay
  - MSPH membrane (Millipore)

**Additional assay technology:** MET Y1230H Lot 001

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



Recombinant Proteins

## ProQinase™ MET Y1230H

Product No.: 0976-0000-1

MET Y1230H Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTLL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQ SMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPEML	KMFDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPOID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHHG	RRRASVAAGI	240
241	<b>LVPRGS</b> PGLD	GICSIEEFKK	<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>AVKSLNRITD</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>VAKGMKYLAS</b>	<b>KKFVHRDLAA</b>	<b>RNCMLDEKFT</b>	<b>VKVADFGLAR</b>	<b>DMHDKKEYYSV</b>	540
541	<b>HNKTGAKLPV</b>	<b>KWMALESLOT</b>	<b>QKFTTKSDVW</b>	<b>SFGVLLWELM</b>	<b>TRGAPPYPDV</b>	<b>NTFDITVYLL</b>	600
600	<b>QGRLLQPEY</b>	<b>CPDPLYEVM</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:** Y1230H

MET wt <sup>1</sup> Amino Acid Sequence							
1	MKAPAVLAPG	ILVLLFTLVQ	RSNGECKEAL	AKSEMNVNMK	YQLPNFTAET	PIQNVILHEH	60
61	HIFLGATNYI	YVLNEEDLQK	VAEYKTGPVL	EHPDCFCQD	CSSKANLSGG	VWKDNINMAL	120
121	VVDYYDDQL	ISCGSVNRGT	QQRHVFPNH	TADIQSEVHC	IFSPQIEEPS	QCPDCVVSAL	180
181	GAKVLSSVKD	RFINFFVGNT	INSSYFPDHP	LHSISVRLK	ETKDGFMFLT	DQSYIDVLP	240
241	FRDSYPIKYV	HAFESNNFIY	FLTVQRETL	AQTFHTRIIR	FCSINSLGHS	YMEMPLECIL	300
301	TEKRKRSTK	KEVFNILQAA	YVSKPGAQLA	RQIGASLNDD	ILFGVFAQSK	PDSAEPMDRS	360
361	AMCAFPKYV	NDFFNKIVNK	NNVRCLQHFY	GNHEHCFNR	TLLRNSSGCE	ARRDEYRTEF	420
421	TTALQRVDF	MGQFSEVLLT	SISTFIKGD	TIANLGTSEG	RFMQVVVSR	GPSTPHVNFL	480
481	LDSPVSPV	IVEHTLNQNG	YTLVITGKI	TKIPLNGLG	RHFQSCSQCL	SAPPFVQCGW	540
541	CHDKCVRSEE	CLSGTWTQOI	CLPAIYKVP	NSAPLEGGTR	LTICGWDFGF	RRNNKFDLKK	600
600	TRVLLGNESC	TLTLESTMN	TLKCTVGPAM	NKHFNMSSII	SNGHGTTQYS	TFSYVDPVIT	660
661	SISPKYGPMA	GGTLLTLTGN	YLNNGNSRHI	SIGGKTCTLK	SVSNSILECY	TPAQTISTEF	720
721	AVKLIKIDLAN	RETSIFSRYE	DPIVYIEHPT	KSFISGGSTI	TGVGKNLNSV	SVPRMVINVH	780
781	EAGRNFTVAC	QHRNSSEIIC	CTTPSLQQLN	LQLPLKTKAF	FMLDGILSKY	FDLIYVHNPV	840
841	FKPFEKPMI	SMGNENVLEI	KGNDIDPEAV	KGEVLKVGNK	SCENIHLHSE	AVLCTVPNDL	900
901	LKLNSELNIE	WKQAISSTVL	GKVIVQPDQN	FTGLIAGVVS	ISTALLLLLG	FFLWLKRRKQ	960
961	<b>IKDLGSELVR</b>	<b>YDARVHTPHL</b>	<b>DRLVSARSVS</b>	<b>PTTEMVSNES</b>	<b>VDYRATFPED</b>	<b>QFPNSSQNGS</b>	1020
1021	<b>CRQVQYPLTD</b>	<b>MSPILTSGDS</b>	<b>DISSPLLQNT</b>	<b>VHIDL SALNP</b>	<b>ELVQAVQHV</b>	<b>IGPSSLIVHF</b>	1080
1081	<b>NEVIGRGHFG</b>	<b>CVYHGTLDN</b>	<b>DGKKIHCAVK</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>GMKYLASKKF</b>	1200
1201	<b>VHRDLAARNC</b>	<b>MLDEKFTVKV</b>	<b>ADFGLARDMY</b>	<b>DKEYYSVH NK</b>	<b>TGAKLPVKWM</b>	<b>ALESLOTQKF</b>	1260
1261	<b>TTKSDVWSFG</b>	<b>VLLWELMTRG</b>	<b>APPYPDVNTF</b>	<b>DITVYLLQGR</b>	<b>RLLQPEYCPD</b>	<b>PLYEVMLK CW</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 fusion protein **Red:** variant in fusion protein

<sup>1</sup>NCBI/Protein accession number NP\_000236.2

**Please notice:**

Variant amino acid counting starting with Ser755 when referring to GenBank accession J02958 (additional 18 aa exon between S755/G756, frequently found in the literature)

This product was manufactured at Reaction Biology in Freiburg, Germany, and is for in vitro research use only, not for use in humans or animals.  
© European Union, 2020. Material may not be reproduced or distributed without written permission from Reaction Biology Europe GmbH