

## DDR2 T654M

discoidin domain receptor tyrosine kinase 2

Recombinant Human Active Protein Kinase

HGNC Symbol: DDR2

Synonyms: MIG20a, NTRKR3, TKT, TYRO10

Product No.: 1193-0000-1

Lot: 001

**Description:** Human DDR2, C-terminal fragment, amino acids R<sub>422</sub>-E<sub>855</sub> (as in [NCBI/Protein](#) entry NP\_006173.2), with a T654M mutation, activated, N-terminal GST-HIS<sub>6</sub> fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

**Product identity:** DDR2 T654M Lot 001, was confirmed as DDR2 by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW<sub>Fusion Protein</sub>:** 77,831 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:** 0.847 µg/µl

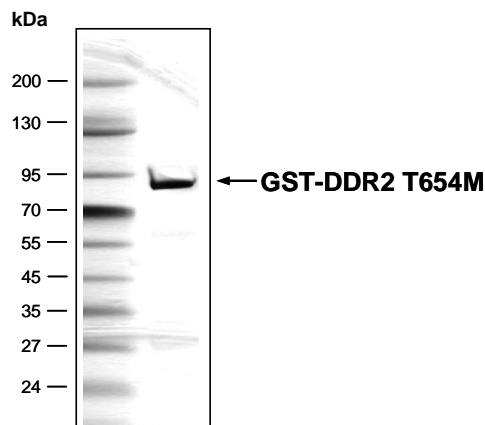
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

### Biochemical Parameters:

Specific kinase activity (P<sub>i</sub> transfer): 13 pmol/µg × min

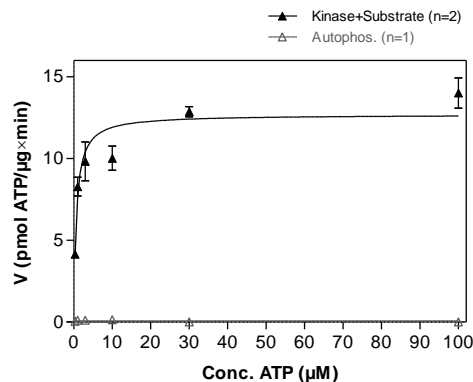
ATP-K<sub>M</sub>: 0.7 µM

### DDR2 T654M Lot 001: Coomassie stain



2.0 µg GST-DDR2 T654M

### DDR2 T654M 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, 80 µg/ml
  - DDR2 T654M: 2 µg/ml
- Filter binding assay
- MSPH membrane (Millipore)

### Additional assay technology:

DDR2 T654M Lot 001 was also successfully tested by ProQinase 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

This product was manufactured at ProQinase in Freiburg, Germany, and is for in vitro research use only, not for use in humans or animals. ProQinase disclaims any warranty explicitly or implied that the use of the product or parts of the product is free from third party intellectual property claims unless this is explicitly stated.

## DDR2 T654M

Product No.: 1193-0000-1

GST-DDR2 T654M Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP EML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI PQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHG	RDSLEVL FQG	240
241	PLAMLRQFWQ	KMLEKASRRM	LDDEMTVSL	LPSDSSMFNN	NRSSSPSEQG	SNSTYDRIFP	300
301	LRPDYQEPSR	LIRKLEFAP	GEEESGCSGV	VKPVQPSGPE	GVPHYAEDI	VNLQGV TGN	360
361	TYSVPAVTMD	LLSGKDVAVE	EFPRKLLTFK	EKLGEQFGE	VHLCEVEGME	KFKDKDFALD	420
421	VSANQPVLVA	VKMLRADANK	NARNDLFKEI	KIMSRLKDPN	IIHLLAVCIT	DDPLCMI MEY	480
481	MENGDLNQFL	SRHEPPNSSS	SDVRTVSYTN	LKFMATQIAS	GMKYLSSLNF	VHRDLATRNC	540
541	LVGKNYTIKI	ADFGMSRNLY	SGDYRIQGR	AVLPIRWSW	ESILLGKFTT	ASDVWAFGVT	600
601	LWETFTFCQE	QPYSQLSDEQ	VIENTGEFFR	DQGRQTYLPQ	PAICPDSVYK	LMLSCWRRDT	660
661	KNRPSFQEIH	LLLLQQGDE					720

1-218: GST Red: HIS6-tag Green: 3C cleavage site blue: DDR2 fragment boxed: T654M mutation

DDR2 wt <sup>1</sup> Amino Acid Sequence							
1	MILIPRMLLV	LFLLLPILSS	AKAQVNPAIC	RYPLGMSGGQ	IPDEDITASS	QWSESTAAY	60
61	GRLDSEEGDG	AWCPEIPVEP	DDLKEFLQID	LHTLHFITLV	GTQGRHAGGH	GIEFAPMYKI	120
121	NYSRDGTRWI	SWRNRHGKQV	LDGNSNPYDI	FLKDLEPIV	ARFVRFIPVT	DHSMNVCMRV	180
181	ELYGCVWLDG	LVSYNAPAGQ	QFVLPGGSI	YLNSVYDGA	VGYSMTEGLG	QLTDGVSGLD	240
241	DFTQTHEYHV	WPGYDYVGR	NESATNGYIE	IMFEFDRIRN	FTTMKVHCNN	MFAKGVKIFK	300
301	EVQCYFRSEA	SEWEPNAISF	PLVLDDVNPS	ARFVTVPLHH	RMASAIKCQY	HFADTWMMFS	360
361	EITFQSDAAM	YNNSEALPTS	PMAPTTYDPM	LKVDDSNTRI	LIGCLVAIIF	ILLAIIVIIL	420
421	WRQFWQKMLE	KASRRMLDDE	MTVSLSLPSD	SSMFNNRNS	SPSEQGSNST	YDRIFPLRPD	480
481	YQEPSRLIRK	LPEFAPGEEE	SGCSGVVKPV	QPSGPEGVPH	YAEADIVNLQ	GVTGGNTYSV	540
541	PAVTMDLLSG	KDVAVEEFPR	KLLTFKEKLG	EGQFGEVHLC	EVEGMEKFKD	KDFALDVSAN	600
601	QPVLVAVKML	RADANKNARN	DFLKEIKIMS	RLKDPNIIHL	LAVCITDDPL	CMITEYMENG	660
661	DLNQFLSRHE	PPNSSSSDVR	TVSYTNLFEM	ATQIASGMKY	LSSLNFVHRD	LATRNC LVGK	720
721	NYTIKIADFG	MSRNLYSGDY	YRIQGRAVLP	IRWMSWESIL	LGKFTTASDV	WAFGVTLWET	780
781	FTFCQE QPYS	QLSDEQVIEN	TGEFFRDQGR	QTYLPQPAIC	PDSVYKMLLS	CWRRDTKNRP	840
841	SFQEIHL LLL	QQGDE					900

blue: DDR2 sequence expressed in recombinant protein Red: variant in recombinant protein

<sup>1</sup>[NCBI/Protein](https://www.ncbi.nlm.nih.gov/Protein/np_006173.2) accession number NP\_006173.2

This product was manufactured at ProQinase in Freiburg, Germany, and is for in vitro research use only, not for use in humans or animals. ProQinase disclaims any warranty explicitly or implied that the use of the product or parts of the product is free from third party intellectual property claims unless this is explicitly stated.