

## ProQinase™ KIT D816V

KIT proto-oncogene receptor tyrosine kinase

Recombinant Human Active Protein Kinase

HGNC Symbol: KIT

Synonyms: CD117, PBT, SCFR, c-Kit

Product No.: 0946-0000-1

Lot: 001

**Description:** Human KIT, C-terminal fragment, amino acids T<sub>544</sub>-V<sub>976</sub> (as in [NCBI/Protein](#) entry NP\_000213.1) with a D816V mutation, N-terminal GST-HIS<sub>6</sub> fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

**Product identity:** KIT D816V Lot 002 was confirmed as KIT by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW**<sub>Fusion Protein</sub>: 77,441 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:** 0.117 µg/µl

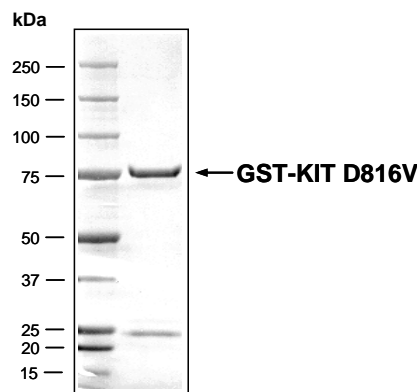
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

**Biochemical Parameters:**

Specific kinase activity (P<sub>i</sub> transfer): 65 pmol/µg × min

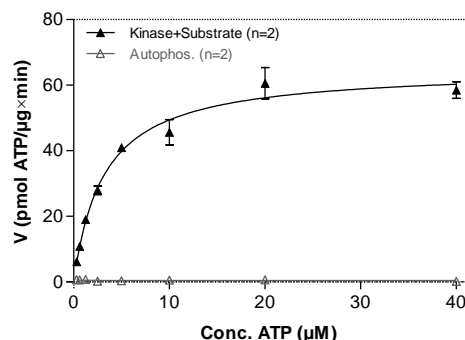
ATP-K<sub>M</sub>: 3.2 µM

**KIT D816V Lot 001:**  
Coomassie stain



2.0 µg GST-KIT D816V

**KIT D816V 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
  - KIT D816V: 4 µg/ml
- Filter binding assay
  - MSPH membrane (Millipore)

**Additional assay technology:**

KIT D816V 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™ KIT D816V

Product No.: 0946-0000-1

| GST-KIT D816V 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  | KRIEAIPOID        | KYLKSSKYIA        | WPLQGWQATF        | GGGDHPPKSD        | PMG <b>HHHHHG</b>  | RDS <b>LEVLFQG</b> | 240 |
| 241  | <b>PLAMGTYKYL</b> | <b>QKPMYEVQWK</b> | <b>VVEEINGNNY</b> | <b>VYIDPTQLPY</b> | <b>DKHWEFPRNR</b>  | <b>LSFGKTLGAG</b>  | 300 |
| 301  | <b>AFGKVVEATA</b> | <b>YGLIKSDAAM</b> | <b>TVAVKMLKPS</b> | <b>AHLTEREALM</b> | <b>SELKVLISYLG</b> | <b>NHMNIVNLLG</b>  | 360 |
| 361  | <b>ACTIGGPTLV</b> | <b>ITEYCCYGD</b>  | <b>LNFLRRKRDS</b> | <b>FICSKQEDHA</b> | <b>EAALYKNLLH</b>  | <b>SKESSCSDST</b>  | 420 |
| 421  | <b>NEYMDMKPGV</b> | <b>SYVVPKADK</b>  | <b>RRSVRIGSYI</b> | <b>ERDVTPAIME</b> | <b>DDELALDLED</b>  | <b>LLSFSYQVAK</b>  | 480 |
| 481  | <b>GMAFLASKNC</b> | <b>IHRDLAARNI</b> | <b>LLTHGRITKI</b> | <b>CDFGLARVIK</b> | <b>NDSNYVVKGN</b>  | <b>ARLPVKWMAE</b>  | 540 |
| 541  | <b>ESIFNCVYTF</b> | <b>ESDVWSYGF</b>  | <b>LWELFSLGSS</b> | <b>PYPGMPVDSK</b> | <b>FYKMIKEGFR</b>  | <b>MLSPEHAPAE</b>  | 600 |
| 601  | <b>MYDIMKTCWD</b> | <b>ADPLKRPTFK</b> | <b>QIVQLIEKQI</b> | <b>SESTNHIYSN</b> | <b>LANCSPNRQK</b>  | <b>PVVDHVSVIN</b>  | 660 |
| 661  | <b>SVGSTASSSQ</b> | <b>PLLVHDDV</b>   |                   |                   |                    |                    | 720 |

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

| KIT wt <sup>1</sup> Amino Acid Sequence |                   |                    |                   |                   |                    |                   |      |
|---|-------------------|--------------------|-------------------|-------------------|--------------------|-------------------|------|
| 1                                       | MARGAWDF          | LCVLLLLLRV         | QTGSSQPSVS        | PGEPSPPSIH        | PGKSDLIVRV         | GDEIRLLCTD        | 60   |
| 61                                      | PGFVKWTFEI        | LDETENENQN         | EWITEKAEAT        | NTGKYTCTNK        | HGLSNSIYVF         | VRDPAKFLV         | 120  |
| 121                                     | DRSLYGKEDN        | DTLVRCPDLD         | PEVTNYSKLG        | CQKPLPKDL         | 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        | AAIAFNVYVN         | TKPEILTYDR        | 420  |
| 421                                     | LVNGMLQCVA        | AGFPEPTIDW         | YFCPGTEQRC        | SASVLPVDVQ        | TLNSSGPPFG         | KLVVQSSIDS        | 480  |
| 481                                     | SAFKHNGTVE        | CKAYNDVGKT         | SAYFNFAFKG        | NNKEQIHPHT        | LFTPLLIGFV         | IVAGMMCIIV        | 540  |
| 541                                     | MILTYKYLQK        | PMYEVQWKVV         | EEINGNNYVY        | IDPTQLPYDH        | KWEFPRNRLS         | FGKTLGAGAF        | 600  |
| 600                                     | <b>GKVVEATAYG</b> | <b>LIKSDAAMTV</b>  | <b>AVKMLKPSAH</b> | <b>LTEREALMSE</b> | <b>LKVLISYLGNH</b> | <b>MNIVNLLGAC</b> | 660  |
| 661                                     | <b>TIGGPTLVIT</b> | <b>EYCCYGDLLN</b>  | <b>FLRRKRDSFI</b> | <b>CSKQEDHAEA</b> | <b>ALYKNLLHSH</b>  | <b>ESSCSDSTNE</b> | 720  |
| 721                                     | <b>YMDMKPGVSY</b> | <b>VVPTKADKRR</b>  | <b>SVRIGSYIER</b> | <b>DVTPAIMEED</b> | <b>ELALDLEDLL</b>  | <b>SFSYQVAKGM</b> | 780  |
| 781                                     | <b>AFLASKNCIH</b> | <b>RDLAARNILL</b>  | <b>THGRITKICD</b> | <b>FGLARDIKND</b> | <b>SNYVVKGNAR</b>  | <b>LPVKWMAPEE</b> | 840  |
| 841                                     | <b>IFNCVYTFES</b> | <b>DVWSYGFIFLW</b> | <b>ELFSLGSSPY</b> | <b>PGMPVDSKPY</b> | <b>KMIKEGFRML</b>  | <b>SPEHAPAEMY</b> | 900  |
| 901                                     | <b>DIMKTCWDAD</b> | <b>PLKRPTFKQI</b>  | <b>VQLIEKQISE</b> | <b>STNHIYSNLA</b> | <b>NCSNPRQKPV</b>  | <b>VDHVSVINRV</b> | 960  |
| 961                                     | <b>GSTASSSQPL</b> | <b>LHVHDDV</b>     |                   |                   |                    |                   | 1020 |

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

<sup>1</sup>[NCBI/Protein](#) accession number NP\_000213.1