

Certificate of Analysis

TGFBR1, active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-912, 14-912-K, 14-912M

Parent Lot # D10EP004N

The data presented in this document apply to the parent lot shown above and to all pack sizes derived from subsequent vialling runs of this parent lot. An alphabetical suffix after the parent lot number is used to denote each vialling run.

Product Description: N-terminal GST-tagged, recombinant, human TGFBR1 amino acids 200–end containing a T204D mutation, expressed by baculovirus in Sf21 insect cells. Purified using glutathione agarose. The T204D mutation has been demonstrated to enhance kinase activity *in vitro* (Wieser *et al*, EMBO (1995);14:2199-2208). Purity 98% by SDS-PAGE and Coomassie blue staining. MW = 62kDa.

Specific Activity (Parent lot# D10EP004N): 273U/mg, where one unit of TGFBR1 activity is defined as 1nmol phosphate incorporated into 2mg/ml casein per minute at 30°C with a final ATP concentration of 100µM.

Formulation: 1.04mg/ml of enzyme in 50mM Tris/HCl pH7.0, 300mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 270mM sucrose, 1mM benzamidine, 0.2mM PMSF, 0.1% 2-mercaptoethanol. Frozen solution.

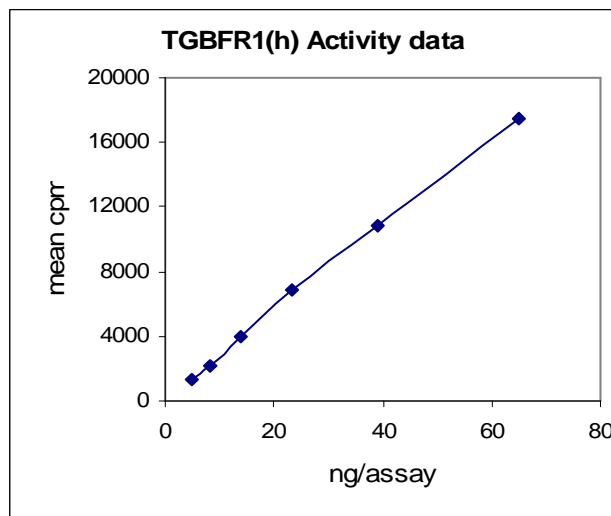
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

Handling Recommendations: Rapidly thaw the vial under cold water and immediately place on ice. Aliquot unused material into pre-chilled microcentrifuge tubes and immediately snap-freeze the vials in liquid nitrogen prior to re-storage at -70°C.

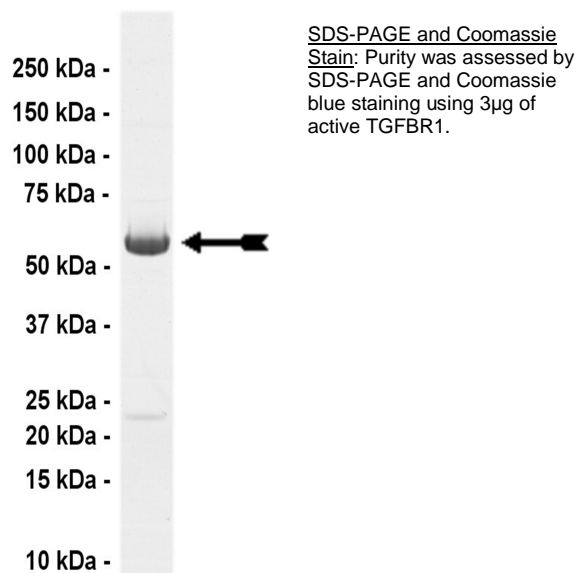
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Kinase Assay: 5–65ng of this lot of enzyme phosphorylated 2mg/ml casein in the assay described on page two. Assay background was subtracted from the actual counts to yield the results shown below.



MS Tryptic Fingerprint: Confirmed identity as TGFBR1 with the translated sequence listed on page three.



Certificate of Analysis

Kinase Assay Protocol

Stock Solutions:

1. **5 x Reaction Buffer:** 40mM MOPS/NaOH pH7.0, 1mM EDTA.
2. **Manganese Chloride:** Use at a final concentration of 1mM. Prepare a 100mM stock and use 0.25µl per well.
3. **TGFBR1, active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 5–65ng per assay point.
4. **[γ -³³P]ATP:** 2.5 x MgAc/[γ ³³P]ATP cocktail: 25mM MgAc and 0.25mM ATP to which is added [γ ³³P]ATP (specific activity approximately 500 - 800cpm/pmol as required).

Assay Procedure (96 well plate format):

1. Add 5µl of 5 x reaction buffer per assay to wells.
2. Add 0.25 µl of 100mM manganese chloride
3. Add 4.75µl of dH₂O.
4. Add 2.5µl of 20mg/ml casein.
5. Add **2.5µl (5–65ng) TGFBR1, active.**
6. Add 10µl of diluted [γ -³³P]ATP mixture.
7. Incubate for 10 minutes at 30°C.
8. Stop the reaction by adding 5µl of 3% phosphoric acid.
9. Transfer a 10µl aliquot onto the appropriate area of a **P30 Filtermat.**
10. Wash the filtermat three times for 5 minutes with 75mM phosphoric acid.
11. Wash the filtermat once for 2 minutes with methanol.
12. Transfer the filtermat to a sealable plastic bag and add 4ml of scintillation cocktail.
13. Read in a scintillation counter. Compare cpm of enzyme samples with cpm of control samples that contain all assay components plus 1µl of 30% phosphoric acid.

Certificate of Analysis

TGFBR1 Sequence Information

<u>Protein</u>	human TGFBR1
<u>Tags</u>	N-terminal GST
<u>Native sequence</u>	T237 of the recombinant protein is equivalent to T200 of human TGFBR1
<u>Accession number</u>	GenBank NM_004612

Recombinant TGFBR1 amino acid sequence:

```

1  MSPILGYWKI  KGLVQPTRL  LEYLEEKYEE  HLYERDEGDK  WRNKKFELGL  EFPNLPYYID
61  GDVKLTQSMA  IIRYIADKHN  MLGGCPKERA  EISMLEGAVL  DIRYGVSRIA  YSKDFETLKV
121  DFLSKLPEML  KMFEDRLCHK  TYLNGDHVTH  PDFMLYDALD  VVLYMDPMCL  DAFPKLVCFK
181  KRIEAIQID   KYLKSSKYIA  WPLQGWQATF  GGGDHPKSD   LEVLFQGPEF  KGLRRQTIAR
241  DIVLQESIGW  GRFGEVWRGK  WRGEEVAVKI  FSSREERSWF  REAEIYQTM   LRHENILGFI
301  AADNKDNGTW  TQLWLVS DYH  EHGSLFDYLN  RYTVTVEGMI  KLALSTASGL  AHLHMEIVGT
361  QGKPAIAHRD  LKSKNILVKK  NGTCCIADLG  LAVRHDSATD  TIDIAPNHRV  GTKRYMAPEV
421  LDD SINMKHF  ESFKRADIYA  MGLVFWEIAR  RCSIGGIHED  YQLPYYDLVP  SDPSVEEMRK
481  VVCEQKLRPN  IPNRWQSCEA  LRVMAKIMRE  CWYANGAARL  TALRIKKTLS  QLSQQEGIKM

```

Recombinant TGFBR1 nucleotide sequence:

```

1  atgtccccta  tactaggtta  ttggaaaatt  aagggccttg  tgcaaccac  tgcacttctt
61  ttggaatata  ttgaagaaaa  atatgaagag  catttgatat  agcgcgatga  aggtgataaa
121  tggcgaaaca  aaaagtttga  attgggtttg  gagtttccca  atcttcctta  ttatattgat
181  ggtgatgtta  aattaacaca  gtctatggcc  atcatacgtt  atatagctga  caagcacaac
241  atgttgggtg  gttgtccaaa  agagcgtgca  gagatttcaa  tgcttgaagg  agcggttttg
301  gatattagat  acgggtgttc  gagaattgca  tatagtaaag  actttgaaac  tctcaaagtt
361  gattttctta  gcaagctacc  tgaaatgctg  aaaatgttct  aagatcgttt  atgtcataaa
421  acatatttaa  atgggatca  tgtaaccat  cctgacttca  tgttgatga  cgctcttgat
481  gttgttttat  acatggacc  aatgtgcctg  gatgcgttcc  caaaattagt  ttgttttaa
541  aaacgtattg  aagctatccc  acaaattgat  aagtacttga  aatccagcaa  gtatatagca
601  tggcctttgc  agggctggca  agccacgttt  ggtggtggcg  accatcctcc  aaaatcggat
661  ctggaagttc  tgttccaggg  gcccgaattc  aaaggcctac  gtcgacaaac  aattgcgaga
721  gatattgtgt  tacaagaaag  cattggcaaa  ggtcgatttg  gagaagtttg  gagaggaaag
781  tggcggggag  aagaagttgc  tgtaagata  ttctcctcta  gagaagaacg  ttcgtggttc
841  cgtgaggcag  agatttatca  aactgtaatg  ttacgtcatg  aaaaatcct  gggatttata
901  gcagcagaca  ataaagaca  tggtaacttg  actcagctct  ggttgggtgc  agattatcat
961  gagcatggat  cccttttga  ttacttaaac  agatacacag  ttactgtgga  agaatgata
1021  aaacttgctc  tgtccacggc  gagcggctct  gcccatcttc  acatggagat  tgttggtagc
1081  caaggaaagc  cagccattgc  tcatagagat  ttgaaatcaa  agaatatctt  ggtaaagaag
1141  aatggaactt  gctgtattgc  agacttagga  ctggcagtaa  gacatgattc  agccacagat
1201  accattgata  ttgctccaaa  ccacagagtg  ggaacaaaaa  ggtacatggc  ccctgaagtt
1261  ctgatgatt  ccataaata  gaaacatttc  gaatccttca  aacgtgctga  catctatgca
1321  atgggcttag  tattctggga  aattgctcga  cgatgttcca  ttggtggaat  tcatgaagat
1381  taccaactgc  cttattatga  tcttgacct  tctgacctat  cagttgaaga  aatgagaaaa
1441  gttgtttgtg  aacagaagtt  aaggccaaat  atcccaaaac  gatggcagag  ctgtgaagcc
1501  ttgagagtaa  tggctaaaat  tatgagagaa  tgttggtagt  ccaatggagc  agctaggctt
1561  acagcattgc  ggattaagaa  aacattatcg  caactcagtc  aacaggaagg  catcaaaatg
1621  taa

```

Reviewed and approved by site quality representative.

Unless otherwise stated in our catalogue or other company documentation accompanying the product(s), our products are intended for research use only and are not to be used for any other purpose, which includes but is not limited to, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses or any type of consumption or application to humans or animals.

© 2014 Eurofins Pharma Discovery Services UK Limited is an independent member of Eurofins Discovery Services.