

Certificate of Analysis

Ret (V804L), active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-758, 14-758-K, 14-758M

Parent Lot # D8CN007U

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 Ret amino acids 658–end containing the V804L mutation. Expressed by baculovirus in Sf21 insect cells. Purified using glutathione agarose.

The V804L mutation has been identified in patients with familial medullary thyroid carcinoma (FMTC) (Lombardo F. et al., (2002), JCEM, 87, 1674-1680; Lesueur F. et al, (2005), JCEM, 90, 3454-3457).

Purity 78% by SDS-PAGE and Coomassie blue staining. MW = 79.2kDa.

Formulation: 1.578mg/ml of enzyme in 50mM Tris/HCl pH7.5, 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.

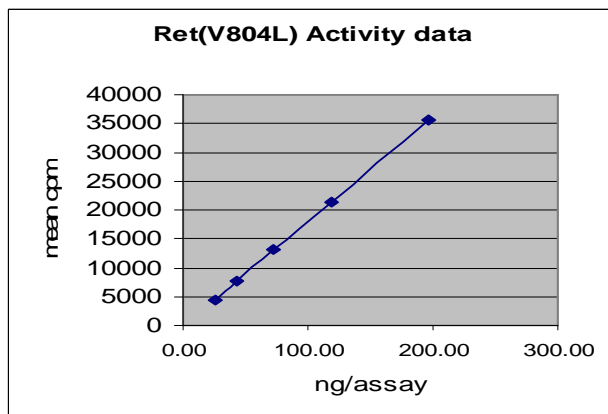
Specific Activity (Parent lot# D8CN007U): 99U/mg, where one unit of Ret(V804L) activity is defined as 1nmol phosphate incorporated into 100µM CHKtide (KKKVSRSGLYRSPSPENLNRPR) per minute at 30°C with a final ATP concentration of 100µM.

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.

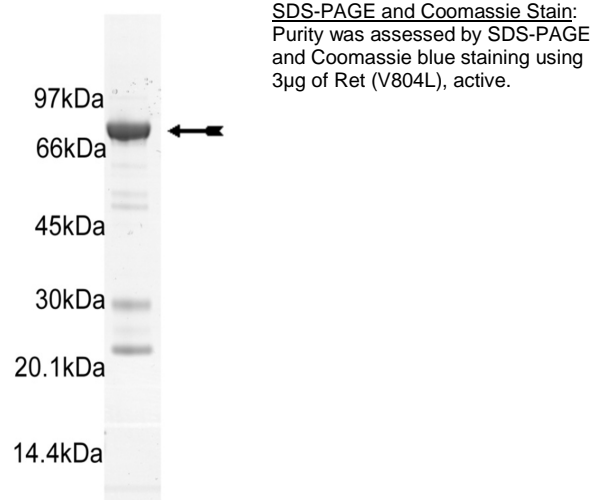
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Kinase Assay: 71–197ng of this lot of enzyme phosphorylated 100µM CHKtide (KKKVSRSGLYRSPSPENLNRPR) 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 Ret(V804L) 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. **CHKtide (KKKVSRSGLYRSPSPENLNRPR):** Use at a final assay concentration of 100 μ M. Prepare a 1mM stock and add 2.5 μ l of stock per assay point.
3. **NaCl:** Use at a final assay concentration of 50mM. Make a 3M stock. Add 0.42 μ l of stock per assay point.
4. **Ret (V804L), active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 5% glycerol, 0.01% Brij-35, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 71–197ng per assay point.
5. **[γ -³³P]ATP:** 2.5 x magnesium acetate/[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 5 μ l of 5 x reaction buffer to wells.
2. Add 2.5 μ l of (KKKVSRSGLYRSPSPENLNRPR).
3. Add 2.5 μ l (71–197ng) Ret(V804L), active.
4. Add 0.42 μ l of 3M NaCl.
5. Add 4.58 μ l of dH₂O.
6. Add 10 μ l of diluted [γ -³³P]ATP mixture.
7. Incubate for 10 minutes at 30°C.
8. Stop the reaction by adding 5 μ l 3% phosphoric acid.
9. Transfer a 10 μ l 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 1 μ l of 30% phosphoric acid.

Certificate of Analysis

Ret (V804L) Sequence Information

<u>Protein</u>	Human Ret (658-end, V804L)
<u>Tags</u>	N-Terminal GST
<u>Native sequence</u>	H237 of recombinant sequence is equivalent to H658 of native human Ret
<u>Accession number</u>	GenBank NM_000323

Recombinant Ret (V804L) amino acid sequence:

```

1 MSPILGYWKI KGLVQPTRLL LEYLEEKYEE HLYERDEGDK WRNKKFELGL EFPNLPYYID
61 GDVKLTQSMA IIRYIADKHN MLGGCPKERA EISMLEGAVL DIRYGVSRIA YSKDFETLKV
121 DFLSKLP EML KMFEDRLCHK TYLNGDHVTH PDFMLYDALD VVLYMDPMCL DAFPKLVCFK
181 KRIEAIQID KYLKSSKYIA WPLQGQWQATF GGGDHPPKSD LEVLFQGPPEF KGLRRQHCHYH
241 KFAHKPPISS AEMTFRRPAQ AFPVSYSSSG ARRPSLDSME NQVSVDAFKI LEDPKWEFPR
301 KNLVLGKTLG EGEFGKVVKA TAFHLKGRAG YTTVAVKMLK ENASPSSELRD LLSEFNVLKQ
361 VNHPHVIKLY GACSQDGPLL LILEYAKYGS LRGFRESRK VGGPYLGS GG SRNSSSLDHP
421 DERALTMGDL ISFAWQISQG MQYLAEMKLV HRDLAARNIL VAEGRKMKIS DFGLSRDVYE
481 EDSYVKRSQG RIPVKWMAIE SLFDHIYTTQ SDVVSFGVLL WEIVTLGGNP YPGIPPERLF
541 NLLKTGHRME RPDNCSEEMY RLMLQCWKQE PDKRPVFADI SKDLEKMMVK RRDYDLAAS
601 TPSDSLIIYDD GLSEETPLV DCNNAPLPRA LPSTWIENKL YGMSDPNWP G ESPVPLTRAD
661 GTNTGFPRYP NDSVYANWML SPSAAKLMDT FDS

```

Recombinant Ret (V804L) nucleotide sequence:

```

1 atgtccccta tactaggtta ttggaaaatt aagggccttg tgcaaccac togacttctt
61 ttggaatata ttgaagaaaa atatgaagag catttgatg agcgcgatga aggtgataaa
121 tggcgaaaca aaaagtttga attgggtttg gagtttcca atcttcctta ttatatgat
181 ggtgatgta aattaacaca gtctatggcc atcatacgtt atatagctga caagcacaac
241 atgttggttg gttgtccaaa agagcgtgca gagatttcaa tgcttgaag agcggttttg
301 gatattagat acgggtgttc gagaattgca tatagtaaag actttgaaac tctcaaagtt
361 gattttctta gcaagctacc tgaaatgctg aaaatgttcg aagatcgttt atgtcataaa
421 acataattaa atggtgatca tgtaacccat cctgacttca tgttgatga cgctcttgat
481 gttgttttat acatggacc aatgtgcctg gatgcttcc caaaatagtt ttgttttaa
541 aaacgtattg aagctatccc acaaattgat aagtacttga aatccagcaa gtatatagca
601 tggcctttgc agggctggca agccacgttt ggtgggtggc accatcctcc aaaatcggat
661 ctggaagttc tgttccaggg gcccgaaatc aaaggcctac gtcgacaaca ctgctaccac
721 aagtttgccc acaagccacc catctcctca gctgagatga ccttccggag gcccgccag
781 gccttccggg tcagctactc ctcttccggt gcccgccggc cctcgctgga ctccatggag
841 aaccaggctc ccgtggatgc cttcaagatc ctggaggatc caaagtggga attcctcgg
901 aagaacttgg ttcttgaaa aactctagga gaaggcgaat ttggaaaagt ggtcaaggca
961 acggccttcc atctgaaagg cagagcaggg tacaccacgg tggcctgtaa gatgctgaaa
1021 gagaacgcct ccccgagtga gcttcgagac ctgctgtcag agttcaacgt cctgaagcag
1081 gtcaaccacc cacatgtcat caaattgtat ggggcctgca gccaggatgg cccgctcctc
1141 ctcatcttgg agtacgcaa atacggctcc ctgctggggt tcctccgca gagccgcaa
1201 gtggggcctg gctacctggg cagtggagcc agccgcaact ccagctcctt ggaccacccg
1261 gatgagcggg ccctcaccat gggcgacctc atctcatttg cctggcagat ctcacagggg
1321 atgcagatc tggccgagat gaagctcggt catcgggact tggcagccag aaacatcctg
1381 gttagctgagg ggcggaagat gaagatttcg gatttcggct tgtcccgaga tgtttatgaa
1441 gaggattcct acgtgaagag gagccagggt cggattccag ttaaattggat ggcaattgaa
1501 tccctttttg atcatatcta caccacgcaa agtgatgat ggtcttttgg tgtcctgctg
1561 tgggagatcg tgaccctagg gggaaacccc tatcctggga ttcctcctga gcggctcttc

```

Certificate of Analysis

```
1621 aaccttctga agaccggcca cgggatggag aggccagaca actgcagcga ggagatgtac
1681 cgctgatgc tgcaatgctg gaagcaggag cggacaaaa ggccggtgtt tgcggacatc
1741 agcaaagacc tggagaagat gatggttaag aggagagact acttggacct tgcggcgtcc
1801 actccatctg actccctgat ttatgacgac ggctctcag aggaggagac accgctgggtg
1861 gactgtaata atgccccctt cctcggagcc ctcccttcca catggattga aaacaaactc
1921 tatggcatgt cagaccgaa ctggcctgga gagagtctg taccactcac gagagctgat
1981 ggcactaaca ctgggtttcc aagatatcca aatgatagtg tatatgctaa ctggatgctt
2041 tcaccctcag cggcaaaatt aatggacacg tttgatagtt aa
```

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.