

## Certificate of Analysis

**CHK2 (I157T), active**  
**(Recombinant enzyme expressed in *E.coli* cells)**  
**Item # 14-741, 14-741-K, 14-741M**  
**Parent Lot # D7DN035U**

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 and C-terminal 6His-tagged recombinant human CHK2, residues 5–end, containing the I157T mutation. Expressed in *E. coli* and purified using glutathione-agarose and Ni<sup>2+</sup>/NTA-agarose.

This I157T mutation is in the forkhead homology-associated (FHA) domain of CHK2 that has been identified in patients with Li-Fraumeni syndrome (a highly penetrant familial cancer phenotype). The I157T mutant retains a similar basal kinase activity compared to the wild type protein. (Wu X *et al*, JBC (2001);276:2971-2974  
 Purity 85.6% by SDS-PAGE and Coomassie staining.  
 MW = 89.6kDa.

**Formulation:** 1.654mg/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.

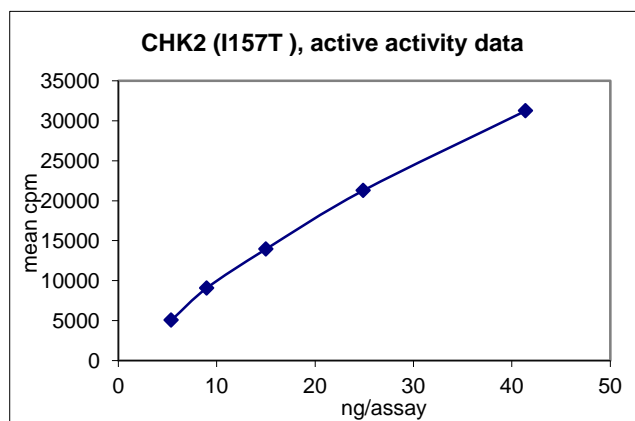
**Specific Activity (Parent lot# D7DN035U):** 867U/mg, where one unit of CHK2(I157T), active activity is defined as 1nmol phosphate incorporated into 100µM CHKtide (KKKVSRSGLYRSPSPENLNRPR) per minute at 30°C with a final ATP concentration of 100µM

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

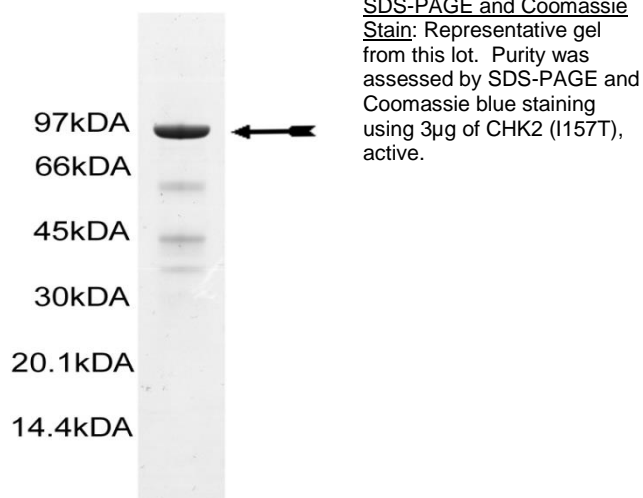
**FOR *IN VITRO* RESEARCH USE ONLY**  
**NOT FOR USE IN HUMANS OR ANIMALS**

### Quality Control Testing

**Kinase Assay:** 3–41ng 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 CHK2, (I157T) with the translated native 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:** Use at a final assay concentration of 100  $\mu$ M. Prepare a 1mM stock and add 2.5 $\mu$ l of stock per assay point.
3. **CHK2, active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 5% glycerol, 0.01% Brij-35, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 3–41ng per assay point.
4. **[ $\gamma$ -<sup>33</sup>P]ATP:** 2.5 x magnesium acetate/[ $\gamma$ -<sup>33</sup>P]ATP cocktail: 25mM MgAc and 0.25mM ATP to which is added [ $\gamma$ -<sup>33</sup>P]ATP (specific activity approximately 500 - 800cpm/pmol as required.)

#### Assay Procedure (96 well plate format):

1. Add 5 $\mu$ l of 5 x reaction buffer per assay to wells.
2. Add 2.5 $\mu$ l of **(KKKVSRSGLYRSPSPENLNRPR)**.
3. Add **2.5 $\mu$ l (3–41ng) CHK2 (I157T) active**.
4. Add 5 $\mu$ l of dH<sub>2</sub>O.
5. Add 10 $\mu$ l of diluted [ $\gamma$ -<sup>33</sup>P] ATP mixture.
6. Incubate for 10 minutes at 30°C.
7. Stop the reaction by adding 5 $\mu$ l of 3% phosphoric acid.
8. Transfer a 10 $\mu$ l aliquot onto the appropriate area of a **P30 Filtermat**.
9. Wash the filtermat three times for 5 minutes with 50mM phosphoric acid.
10. Wash the filtermat once for 2 minutes with methanol.
11. Transfer the filtermat to a sealable plastic bag and add 4ml of scintillation cocktail.
12. Read in a scintillation counter. Compare cpm of enzyme samples with cpm of control samples that contain all assay components plus 1 $\mu$ l of 30% phosphoric acid.

## Certificate of Analysis

### CHK2 (I157T) Sequence Information

<b><u>Protein</u></b>	Human CHK2 (5-end, I157T)
<b><u>Tags</u></b>	N-terminal GST and C-terminal 6His tags
<b><u>Native sequence</u></b>	S249 of recombinant sequence is equivalent to S5 of native human CHK2
<b><u>Accession number</u></b>	GenBank NM 007194

#### Recombinant CHK2 amino acid sequence:

```

1 MSPISRMPIL GYWKIKGLVQ PTRLLLEYLE EKYEEHLYER DEGDKWRNKK FELGLEFPNL
61 PYYIDGDVKL TQSMAIIRYI ADKHNMLGGC PKERAEISML EGAVLDIRYG VSRIAYSKDF
121 ETLKVDFLSK LPEMLKMFED RLCHKTYLNG DHVTHPDFML YDALDVVLYM DPMCLDAFPK
181 LVCFKKRIEA IPQIDKYLKS SKYIAWPLQG WQATFGGGDH PPKSDLVPRG SRRASVGSJM
241 PMSRPRRPSD VEAQQSHGSS ACSQPHGSVT QSQSSSSQSQ GISSSSTSTM PNSSQSSHSS
301 SGTLSSETV STQELYSIPE DQEPEDQEPE EPTPAPWARL WALQDGFANL ECVNDNYWFG
361 RDKSCEYCFD EPLLKRTDKY RTYSKKHFRI FREVGPKNSY TAYIEDHSGN GTFVNTLVG
421 KGKRRPLNNN SEIALSLSRN KVFVFFDLTV DDQSVYPKAL RDEYIMSKTL GSGACGEVKL
481 AFERKTCKKV AIKIISKRFK AIGSAREADP ALNVETEIEI LKKNLHPCII KIKNFFDAED
541 YYIVLELMEG GELFDKVVGN KRLKEATCKL YFYQMLLAVQ YLHENGIIHR DLKPENVLLS
601 SQEEDCLIKI TDFGHSKILG ETSLMRTL CG TPTYLAPEVL VSVGTAGYNR AVDCWSLGI
661 LFICLSGYPP FSEHRTQVSL KDQITSGKYN FIPEVWAEVS EKALDLVKKL LVVDPKARFT
721 TEEALRHPWL QEDMKRKFQ DLLSEENEST ALPQVLAQPS TSRKRPREGE AEGAETTKRP
781 AVCAAVLHHH HHH

```

#### Recombinant CHK2 nucleotide sequence:

```

1 atgtccccta tatctagaat gcctatacta gggttattgga aaattaaggg ccttgtgcaa
61 cccactcgac ttccttttga atactctgaa gaaaaatatg aagagcattt gtatgagcgc
121 gatgaagggtg ataaatggcg aaacaaaaag tttgaattgg gtttggagtt tcccaatctt
181 ccttattata ttgatggatga tgttaaatta acacagtcta tggccatcat acgttatata
241 gctgacaagc acaacatggt gggtgggtgt ccaaaagagc gtgcagagat ttcaatgctt
301 gaaggagcgg ttttggatat tagatacggg gtttcgagaa ttgcatatag taaagacttt
361 gaaactctca aagttgattt tcttagcaag ctacctgaaa tgctgaaaaa gttcgaagat
421 cgtttatgct ataaaacata tttaaatggt gatcatgtaa cccatcctga cttcatgttg
481 tatgacgctc ttgatgttgt tttatacatg gacccaatgt gcctggatgc gttcccaaaa
541 ttagtttggg ttataaaacg tattgaagct atcccacaaa ttgataagta cttgaaatcc
601 agcaagtata tagcatggcc tttgcaagggc tggcaagcca cgtttgggtg tggcgacct
661 ctcccaaaat cggatctggt tccgcgtgga tctcgtcgtg catctgttgg atcgcatatg
721 cccatgtcga ggccacgaag gccgtcggat gttgaggctc agcagtctca tggcagcagt
781 gcctgttcac agccccatgg cagcgttacc cagtcccaag gctcctcctc acagtcccag
841 ggcatatcca gctcctctac cagcacgatg ccaaactcca gccagtctc tcaactccagc
901 tctgggacac tgagctcctt agagacagtg tccactcagg aactctattc tattcctgag
961 gaccaagaac ctgaggacca agaactgag gaggcctacc ctgccccctg ggctcgatta
1021 tgggcccttc aggatggatt tgccaatctt gaatgtgtga atgacaacta ctggtttggg
1081 agggacaaaa gctgtgaata ttgctttgat gaaccactgc tgaagaagac agataaatac
1141 cgaacataca gcaagaaaca ctttcggatt ttcagggaag tgggtcctaa aaactcttac
1201 actgcataca tagaagatca cagtggcaat ggaacctttg taaatacaga gctttagtagg
1261 aaaggaaaac gccgtccttt gaataacaat tctgaaattg cactgtcact aagcagaaat
1321 aaagtttttg tcttttttga tctgactgta gatgatcagt cagtttatcc taaggcatta
1381 agagatgaat acatcatgct aaaaactctt ggaagtgggt cctgtggaga ggtaaagctg
1441 gctttcgaga ggaaaacatg taagaaagta gccataaaga tcatcagcaa aaggaagttt

```

## Certificate of Analysis

```
1501 gctattgggtt cagcaagaga ggcagaccga gctctcaatg ttgaaacaga aatagaaatt
1561 ttgaaaaagc taaatcatcc ttgcatcatc aagattaata acttttttga tgcagaagat
1621 tattatattg ttttgggaatt gatggaaggg ggagagctgt ttgacaaagt ggtggggaat
1681 aaacgcctga aagaagctac ctgcaagctc ttttttacc agatgctctt ggctgtgcag
1741 taccttcatg aaaacggtat tatacaccgt gacttaaagc cagagaatgt tttactgtca
1801 tctcaagaag aggactgtct tataaagatt actgattttg ggcactcaa gattttggga
1861 gagacctctc tcatgagaac cttatgtgga acccccacct acttggcgcc tgaagtctt
1921 gtttctgttg ggactgctgg gtataaccgt gctgtggact gctggagttt aggagttatt
1981 ctttttatct gccttagtgg gtatccacct ttctctgagc ataggactca agtgtcactg
2041 aaggatcaga tcaccagtgg aaaatacaac ttcattcctg aagtctgggc agaagtctca
2101 gagaaagctc tggaccttgt caagaagttg ttggtagtgg atccaaaggc acgttttacg
2161 acagaagaag ccttaagaca cccgtggctt caggatgaag acatgaagag aaagtttcaa
2221 gatcttctgt ctgaggaaaa tgaatccaca gctctacccc aggttctagc ccagccttct
2281 actagtcgaa agcggccccg tgaaggggaa gccgagggtg ccgagaccac aaagcgccca
2341 gctgtgtgtg ctgctgtgtt gcatcaccat caccatcact ga
```

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