

Certificate of Analysis

PKB beta (S474D)

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-339, 14-339-K, 14-339M

Parent Lot # D8JN055U

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 6His-tagged recombinant human PKB beta residues 120–end containing the S474D mutation. Expressed by baculovirus in Sf21 insect cells. Purified using Ni²⁺/NTA agarose, activated by PDK1, and repurified by heparin-Sepharose. Purity 81.1% by SDS-PAGE and Coomassie blue staining. MW = 42.8kDa.

Specific Activity (Parent lot# D8JN055U): 276U/mg, where one unit of PKB beta activity is defined as 1nmol phosphate incorporated into 30µM modified crosstide (GRPRTSSFAEGKK) per minute at 30°C with a final ATP concentration of 100µM.

Formulation: 0.662mg/ml of enzyme in 50mM Tris/HCl pH7.5, 270mM sucrose, 150mM NaCl, 1mM EGTA, 1mM benzamidine, 0.2mM PMSF, 0.1 % 2-mercaptoethanol, 0.03 % Brij-35. Frozen solution.

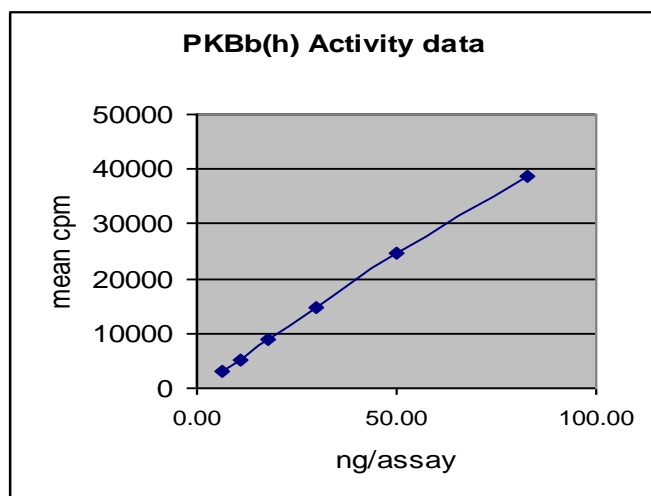
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

Handling Recommendations: Rapidly thaw the vial under cold water and immediately place on ice. Aliquot unused material into pre-chilled 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: 6.6–82.8ng of this lot of enzyme phosphorylated 30µM modified crosstide 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 PKB beta 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. **Modified Crosstide (GRPTSSFAEGKK):** Use at a final assay concentration of 30 μ M. Prepare a 300 μ M stock and add 2.5 μ l of stock per assay point.
3. **PKB beta, active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 5% glycerol, 0.01% Brij-35, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 6.6–82.8ng per assay point.
4. **[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 per well.
2. Add 2.5 μ l of **modified crosstide**.
3. Add **2.5 μ l (6.6–82.8ng) PKB beta, active**.
4. Add 5 μ l of dH₂O.
5. Add 10 μ l of diluted [γ -³³P]ATP mixture.
6. Incubate for 10 minutes at 30°C.
7. Stop the reaction by adding 5 μ l of 3% phosphoric acid.
8. Transfer a 10 μ 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 μ l of 30% phosphoric acid.

Certificate of Analysis

PKB beta Sequence Information

Protein	human PKB beta (S474D)
Tags	N-terminal 6His
Native sequence	M8 of the recombinant protein is equivalent to M120 of human PKB beta
Accession number	EMBL M95936. The recombinant protein contains an amino acid substitution, S474D, with reference to the native sequence. This mimics the phosphorylation of S474.

Recombinant PKB beta amino acid sequence:

```

1  MHHHHHHMDY KCGSPSDSST TEEMEVAVSK ARAKVTMNDF DYLLKLLGKGT FGKVILVREK
61  ATGRYYAMKI LRKEVIIAKD EVAHTVTESR VLQNTRHPFL TALKYAFQTH DRLCFVMEYA
121 NGGELFFHLS RERVFTTEERA RFYGAEIVSA LEYLHSRDVV YRDIKLENLM LDKDGHKIT
181 DFGLCKEGIS DGATMKTFCTG TPEYLAPEVL EDNDYGRAVD WWGLGVVMEY MMCGRLPFYN
241 QDHERLFELI LMEEIRFPRT LSPEAKSLLA GLLKKDPKQR LGGGPSDAKE VMEHRFFLSI
301 NWQDVVQKKL LPPFKPQVTS EVDTRYFDDE FTAQSITITP PDRYDSLGLL ELDQRTTHFPQ
361 FDYSASIRE

```

Recombinant PKB beta nucleotide sequence:

```

1  atgcatcatc accatcacca tatggactac aagtgtggct cccccagtga ctctccacg
61  actgaggaga tggaagtggc ggtcagcaag gcacgggcta aagtgaccat gaatgacttc
121 gactatctca aactccttgg caagggaacc tttggcaaag tcatcctggt gcgggagaag
181 gccactggcc gctactacgc catgaagatc ctgcggaagg aagtcatcat tgccaaggat
241 gaagtcgctc acacagtcac cgagagccgg gtccctccaga acaccaggca cccgttcctc
301 actgcgctga agtatgcctt ccagaccacac gaccgcctgt gctttgtgat ggagtatgcc
361 aacgggggtg agctgttctt ccacctgtcc cgggagcgtg tcttcacaga ggagcggggc
421 cggttttatg gtgcagagat tgtctcggct cttgagtact tgcactcgcg ggacgtggta
481 taccgcgaca tcaagctgga aaacctcatg ctggacaaag atggccacat caagatcact
541 gactttggcc tctgcaaaga gggcatcagt gacggggcca ccatgaaaac cttctgtggg
601 accccggagt acctggcgcc tgaggtgctg gaggacaatg actatggccg ggccgtggac
661 tgggtggggc tgggtgtggt catgtacgag atgatgtgcg gccgcctgcc cttctacaac
721 caggaccacg agcgctctt cgagctcatc ctcatggaag agatccgctt cccgcgcacg
781 ctgagccccg aggccaaagtc cctgcttgct gggctgctta agaaggacct caagcagagg
841 cttggtgggg ggcccagcga tgccaaggag gtcattggagc acaggttctt cctcagcatc
901 aactggcagg acgtggtcca gaagaagctc ctgccaccct tcaaacctca ggtcacgtcc
961 gaggtcgaca caaggtactt cgatgatgaa tttaccgccc agtccatcac aatcacacct
1021 cctgaccgct atgacagcct gggcttactg gagctggacc agcggaccca cttccccag
1081 ttcgactact cggccagcat ccgcgagtga

```

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.