

Certificate of Analysis

ULK1, active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-959, 14-959-K, 14-959M

Parent Lot # D15BP002N

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 ULK1 amino acids 1-314 expressed by baculovirus in Sf21 insect cells. Purified using immobilized metal affinity chromatography followed by gel filtration. Purity 83% by SDS-PAGE and Coomassie blue staining. MW = 39kDa.

Specific Activity (Parent lot# D15BP002N): 2858U/mg, where one unit of ULK1 activity is defined as 1nmol phosphate incorporated into 2mg/ml casein per minute at 30°C with a final ATP concentration of 100µM.

Formulation: 0.45mg/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.

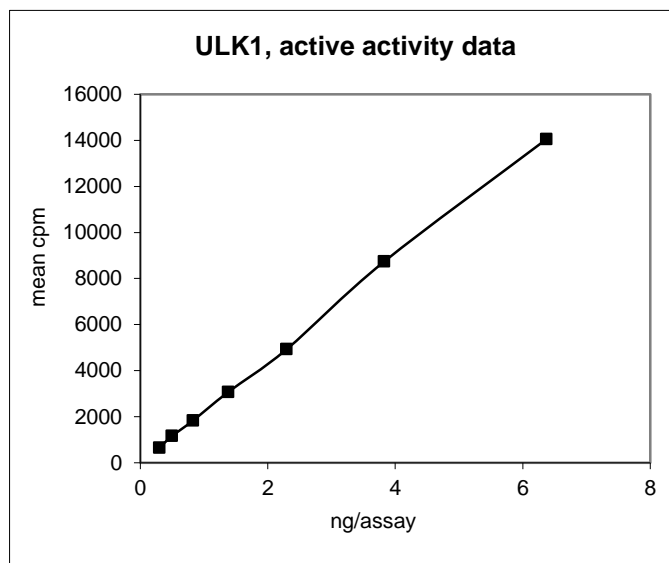
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

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.

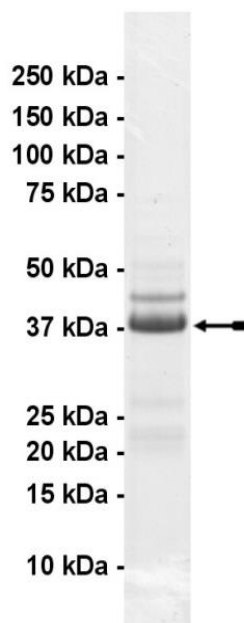
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Kinase Assay: 0.3–6.4ng 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 ULK1 with the translated sequence listed on page three



SDS-PAGE and Coomassie Stain:
Purity was assessed by SDS-PAGE and Coomassie blue staining using 3µg of ULK1, active

Certificate of Analysis

Kinase Assay Protocol

Stock Solutions:

1. **5 x Reaction Buffer:** 40mM MOPS/NaOH pH7.0, 1mM EDTA.
2. **Casein:** Use at a final assay concentration of 2mg/ml. Prepare a 20mg/ml stock and add 2.5µl of stock per assay point.
3. **ULK1, active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 0.3–6.4ng 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 2.5µl of 2mg/ml casein.
3. Add **2.5µl (0.3–6.4ng) ULK1, active.**
4. Add 5µl of dH₂O.
5. Add 10µl of diluted [γ -³³P]ATP mixture.
6. Incubate for 10 minutes at 30°C.
7. Stop the reaction by adding 5µl of 3% phosphoric acid.
8. Transfer a 10µl aliquot onto the appropriate area of a **P30 Filtermat.**
9. Wash the filtermat three times for 5 minutes with 75mM 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 1µl of 30% phosphoric acid.

Certificate of Analysis

ULK1, active Sequence Information

Protein	human ULK1
Tags	N-terminal 6His
Native sequence	M31 of the recombinant protein is equivalent to M1 of human ULK1.
Accession number	GenBank NM_003565.1

Recombinant ULK1 amino acid sequence:

```

1 MSYYHHHHHH DYDIPTTENL YFQGAMDPEF MEPGRGGTET VGKFEFSRKD LIGHGAFVAV
61 FKGRHREKHD LEVAVKCINK KNLAKSQTLG GKEIKILKEL KHENIVALYD FQEMANSVYL
121 VMEYCNGGDL ADYLHAMRTL SEDTIRLFLQ QIAGAMRLLH SKGIIHRDLK PQNILLSNPA
181 GRRANPNSIR VKIADFGFAR YLQSNMMAAT LCGSPMYMAP EVIMSQHYDG KADLWSIGTI
241 VYQCLTGKAP FQASSPQDLR LFYEKNKTLV PTIPRETSAP LRQLLLALLQ RNHKDRMDFD
301 EFFHHPFLLD SPSVRKSPPV PVPSYPSSGS GSSSSSSSTS HLAS

```

Recombinant ULK1 nucleotide sequence:

```

1 atgtcgtact accatcacca tcaccatcac gattacgata tcccaacgac cgaaaacctg
61 tatttttcagg gcgccatgga tccggaattc atggagcccg gccgcggcgg cacagagacc
121 gtgggcaagt tcgagttctc ccgcaaggac ctgatcggcc acggcgcctt cgcggtggtc
181 ttcaagggcc gccaccgcca gaagcacgat ttggaggctc ccgtcaagtg cattaacaag
241 aagaacctcg ccaagtctca gacgtgctg gggaaggaaa tcaaaatcct gaaggaactg
301 aaacatgaaa acatcgtggc cctgtacgac ttccaggaaa tggctaattc tgtctacctg
361 gttatggagt actgcaacgg tggggacctg gccgactacc tgcacgcat gcgcacgctg
421 agcgaggaca ccacaggct cttcctgcag cagatcgcg gccccatgcg gcttctgcac
481 agcaaaggca tcatccaccg cgacctgaaa ccgcagaaca tcctgctgtc caaccccgcc
541 ggccgcccgc ccaaccccaa cagcatccgc gtcaagatcg ctgacttcgg cttcgcgcg
601 tacctccaga gcaacatgat ggcggccaca ctctgcggct ccccatgta catggcccc
661 gaggtcatca tgtcccagca ctacgacggg aaggcggacc tgtggagcat cggcaccatc
721 gtctaccagt gcctgacggg gaaggcgccc ttccaggcca gcagcccca ggacctgcgc
781 ctgttctacg agaagaacaa gacgttggtc cccaccatcc cccgggagac ctcgcccccg
841 ctgcggcagc tgctcctggc cctactgcaa cgcaaccaca aggaccgat ggacttcgat
901 gagttttttc atcacccttt cctcgatgcc agcccctcgg tcaggaaatc cccaccctg
961 cctgtgcct cgtaccaag ctcggggtcc ggcagcagct ccagcagcag ctccacctcc
1021 cacctggcct cctaa

```

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