

LANCE® Ultra

**ULight™ - labeled Acetyl-CoA carboxylase (Ser79) Peptide****Product number:** TRF0133-M **Lot Number:** 3221372**Product Format:** TRF0133-D: 0.5 nmole (100µL)  
TRF0133-M: 5 nmoles (Lyophilized, add 1mL of water)**Manufacturing date:** 11/20/2023 **Document version:** 1**Product Information**

**Phosphorylation Motif:** RSAM<sub>S</sub>GLHL  
Synthetic peptide derived from residues 73-85 of rat Acetyl CoA Carboxylase (also known as SAMS peptide) in which Ser77 is mutated to Ala; phosphorylation site: Ser79

**Storage Buffer:** 50 mM Tris-HCl (pH 7.4), 0.9% NaCl, 0.1% BSA and 0.05% sodium azide as preservative

**Molecular Weight:** 2 674

**Stability:** This product is stable for at least **9 months** from the manufacturing date when stored in its original packaging and the recommended storage conditions.

**Storage Conditions:** Store at -20°C. Store protected from light. Repeated freezing and thawing should be avoided.

**Quality Control**

The QC release specifications are based on spectrophotometric analysis of the labeled protein. We certify that results meet our quality release criteria.

**Labeling Ratio:** 1.0 (dye molecule/peptide)  
**Concentration:** 13.4 µg/mL (5 µM)

**Recommended Assay Conditions****AMPK $\alpha$ 1 kinase: ATP titration**

## SUGGESTED METHOD:

*(Specific applications might require optimization)*Reagent Preparation

- Prepare 1X Kinase Assay Buffer: 50 mM HEPES pH 7.5, 1 mM EGTA, 10 mM MgCl<sub>2</sub>, 2 mM DTT and 0.01% Tween-20.
- Prepare a 2X AMPK $\alpha$ 1 solution: dilute enzyme to a concentration of 4 nM in Kinase Assay Buffer. Keep on ice.
- Prepare a 4X ULight-Acetyl CoA Carboxylase (Ser79) peptide solution: dilute ULight-Acetyl CoA Carboxylase (Ser79) to a concentration of 200 nM in Kinase Assay Buffer.
- Prepare a 4X ATP solution: dilute ATP to concentrations ranging from 40 nM to 4 mM (serial half-log dilutions) in Kinase Assay Buffer. Keep on ice.
- Prepare 1X Detection Buffer: dilute 0.5 mL of 10X Detection Buffer with 4.5 mL of H<sub>2</sub>O.
- Prepare a 4X Stop solution\*: dilute EDTA to a concentration of 40 mM in 1X Detection Buffer.
- Prepare a 4X Detection Mix: dilute Europium-anti-phospho-Acetyl CoA Carboxylase (Ser79) antibody to a concentration of 8 nM in 1X Detection Buffer.

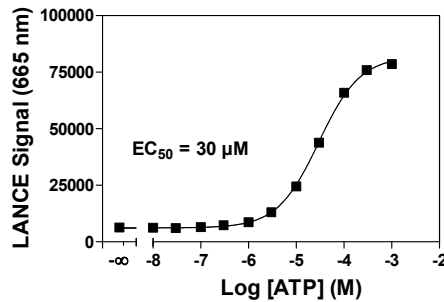
Protocol

- Pipet 5  $\mu$ L of 2X AMPK $\alpha$ 1 solution into a 384-well white OptiPlate-384 (2 nM final concentration).
- Add 2.5  $\mu$ L of 4X ULight- Acetyl CoA Carboxylase (Ser79) solution (50 nM final concentration).
- Add 2.5  $\mu$ L of 4X ATP solution (10 nM to 1 mM final concentrations).
- Cover plate with TopSeal-A and incubate 30 min at 23°C.
- Add 5  $\mu$ L of 4X Stop solution\* and incubate 5 min at 23°C.
- Add 5  $\mu$ L of 4X Detection Mix (2 nM Eu-anti-phospho-Acetyl CoA Carboxylase (Ser79) final conc.).
- Cover plate with TopSeal-A and incubate 60 min at 23°C.
- Remove TopSeal-A and read in TR-FRET mode at 665 nm (excitation at 320 or 340 nm).

\*Alternatively, the Stop solution and Detection Mix can be premixed and added together to the kinase reaction.

## Typical Product Data

AMPK $\alpha$ 1 kinase assay using ULight-Acetyl CoA Carboxylase (Ser79) Peptide and Eu-anti-phospho-Acetyl CoA Carboxylase (Ser79) Antibody obtained using the EnVision® Multilabel Reader:



## Suggested Materials

	Supplier
• Substrate: ULight™ - Acetyl CoA Carboxylase (Ser79)Pept.	Revvity Inc
• Antibody: Eu- anti- phospho-Acetyl CoA carboxylase(Ser79)	Revvity Inc
• Kinase: AMPK $\alpha$ 1	Carna Biosciences
• Detection Buffer: LANCE® Detection Buffer, 10X	Revvity Inc
• Plate: OptiPlate™-384, white	Revvity Inc
• TopSeal™-A	Revvity Inc

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

TRF0133-R Rev01

revvity

Revvity, Inc.  
940 Winter Street  
Waltham, MA 02451 USA

(800) 762-4000 [www.revvity.com](http://www.revvity.com)

For a complete listing of our global offices, visit [www.revvity.com](http://www.revvity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.