

LANCE® Ultra

Eu-W1024 labeled Anti-phospho Acetyl-CoA carboxylase (Ser79)**Antibody**

Product number: TRF0208-D **Lot Number:** 3254971

Product Format: TRF0208-D: 10 µg
TRF0208-M: 100 µg

Manufacturing date: 8/23/2023 **Document version:** 1

Product Information

Antibody: Europium-labeled mouse monoclonal antibody recognizing phospho-Ser79 in rat acetyl CoA carboxylase.

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

Molecular Weight: 160 000

Stability: This product is stable for at least 24 months from the manufacturing date when stored in its original packaging and at the recommended storage conditions.

Storage Conditions: Store at 4°C

Quality Control

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

Labeling Ratio: 7.5
Concentration: 100 µg/mL (0.625 µM)

Recommended Assay Conditions**AMPK α 1 KINASE: ATP TITRATION**

SUGGESTED METHOD:

(Specific applications might require optimization)

Reagent Preparation

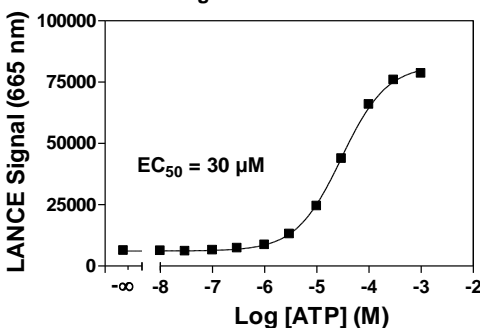
- Prepare 1X Kinase Assay Buffer: 50 mM HEPES pH 7.5, 1 mM EGTA, 10 mM MgCl₂, 2 mM DTT and 0.01% Tween-20.
- Prepare a 2X AMPK α 1 solution: dilute enzyme to a concentration of 4 nM in Kinase Assay Buffer. Keep on ice.
- Prepare a 4X ULight-Acetyl CoA Carboxylase (Ser79) peptide solution: dilute ULight-Acetyl 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 H₂O.
- 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 μ L of 2X AMPK α 1 solution into a 384-well white OptiPlate-384 (2 nM final concentration).
 - Add 2.5 μ L of 4X ULight-Acetyl CoA carboxylase (Ser79) solution (50 nM final concentration).
 - Add 2.5 μ 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 μ L of 4X Stop solution* and incubate 5 min at 23°C.
 - Add 5 μ 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 α 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

- | | |
|---|-------------------|
| • Substrate: ULight [™] - Acetyl CoA Carboxylase (Ser79) Pept. | Supplier |
| • Antibody: Eu- anti-phospho-Acetyl CoA Carboxylase (Ser79) | Revvity Inc |
| • Kinase: AMPK α 1 | Revvity Inc |
| • Detection Buffer: LANCE [®] Detection Buffer, 10X | Carna Biosciences |
| • Plate: OptiPlate [™] -384, white | Revvity Inc |
| | 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.

TRF0208-R Rev01

The logo for Revvity, featuring the word "revvity" in a lowercase, sans-serif font.

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

(800) 762-4000 www.revvity.com

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2023, Revvity, Inc. All rights reserved.