

AMPK α 1 Kinase Assay

ULight-Acetyl-CoA Carboxylase (Ser79) Peptide and Europium-anti-phospho-Acetyl-CoA Carboxylase (Ser79) Antibody

Two LANCE Ultra companion products – two convenient sizes!

ULight™-Acetyl-CoA Carboxylase (Ser79) Peptide (SAMS Peptide):

- TRF0118-D: 0.5 nmole, 1,000* assay points
- TRF0118-M: 5 nmoles, 10,000* assay points

*0.5 pmol/assay point

PEPTIDE SEQUENCE:

CHMRSAMSSGLHLVKRR

Synthetic peptide derived from residues 73-85 of rat acetyl-CoA carboxylase in which Ser77 is mutated to Ala; phosphorylation site: Ser79.

VALIDATED FOR KINASE: AMPK α 1

POTENTIAL SUBSTRATE FOR KINASES:

AMP-activated subfamily of protein kinases

Europium-anti-phospho-Acetyl-CoA Carboxylase (Ser79) Antibody:

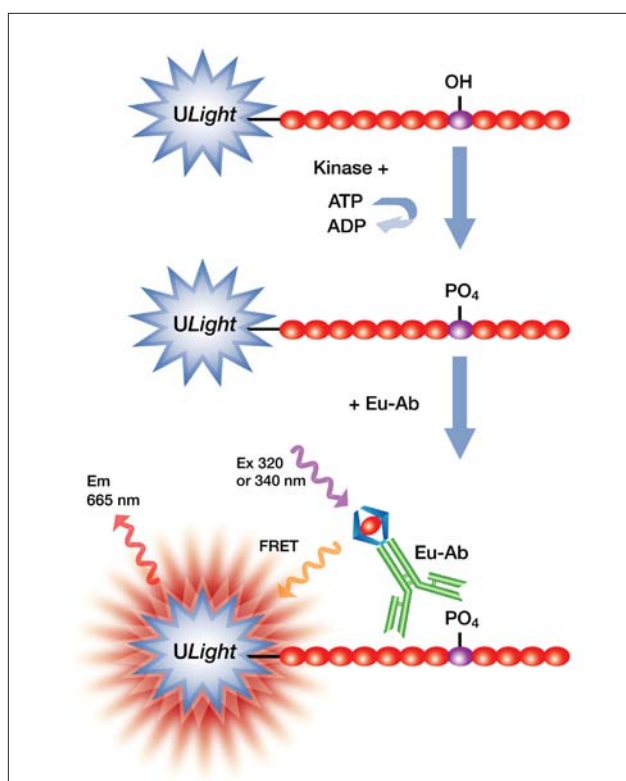
- TRF0208-D: 10 μ g, 1,562* assay points
- TRF0208-M: 100 μ g, 15,625* assay points

*40 fmol/assay point

RECOGNIZED MOTIF:

SSMpSGL

Europium-labeled mouse monoclonal antibody recognizing phospho-Ser79 of rat acetyl-CoA carboxylase.



LANCE Ultra Kinase Assays

LANCE® Ultra time-resolved fluorescence resonance energy transfer (TR-FRET) assays use a proprietary europium chelate donor dye, W-1024 (Eu), with ULight, an innovative small molecular weight acceptor dye with a red-shifted fluorescent emission. In kinase assays, the binding of a Eu-labeled anti-phospho-substrate antibody to the phosphorylated ULight-substrate brings donor and acceptor molecules into close proximity.

After irradiation of the kinase reaction at 320 or 340 nm, the energy from the Eu donor is transferred to the ULight acceptor dye which, in turn, generates light at 665 nm. The intensity of the light emission is proportional to the level of ULight-substrate phosphorylation.

Development of an AMPK α 1 Kinase Assay

Additional reagents

AMPK α 1 active	Carna # 02-113
LANCE Detection Buffer, 10X	PerkinElmer # CR97-100
OptiPlate™-384, white	PerkinElmer # 6007299
TopSeal™-A	PerkinElmer # 6005185
Kinase Buffer: 50 mM HEPES pH 7.5, 1 mM EGTA, 10 mM MgCl ₂ , 2 mM DTT and 0.01% Tween-20.	

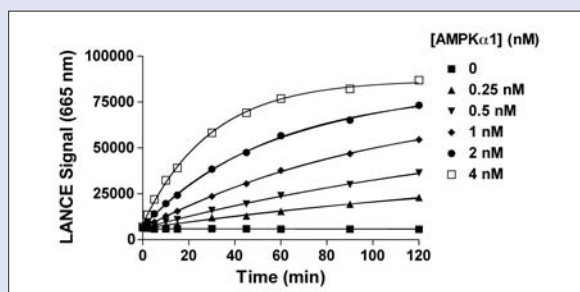
Suggested procedure

- Dilute the AMPK α 1 kinase, ATP, inhibitors and *ULight*-Acetyl-CoA Carboxylase peptide in Kinase Buffer.
- Prepare a 4X Detection Mix by diluting the Eu-anti-phospho Acetyl-CoA Carboxylase antibody to 8 nM in 1X LANCE Detection Buffer.
- Add to the wells of a white Optiplate-384:
 - 5 μ L of AMPK α 1 enzyme
 - 2.5 μ L of inhibitor or Kinase Buffer
 - 2.5 μ L of *ULight*-Acetyl-CoA Carboxylase peptide/ATP mix (for ATP titration, ATP dilutions are added separately in Kinase Buffer).

- Cover the plate with TopSeal-A and incubate at room temperature (RT).
- Stop kinase reactions by adding 5 μ L of 40 mM EDTA prepared in 1X Detection Buffer (Stop Solution). Leave for 5 min at RT.
- Add 5 μ L of 4X Detection Mix (Eu-anti-phospho-Acetyl-CoA Carboxylase Antibody at a final concentration of 2 nM).
- Cover with TopSeal-A and incubate for 1 h at RT.
- Remove TopSeal-A and read signal with the EnVision[®] Multilabel Reader in TR-FRET mode (excitation at 320 nm and emission at 665 nm).

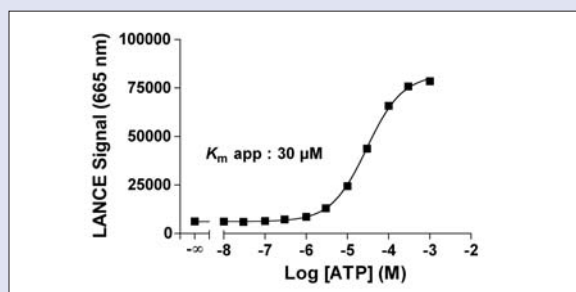
NOTE: Eu-labeled antibodies and EDTA can be premixed before use as a 2X concentrated Stop Solution/Detection Mix to minimize the number of liquid handling steps.

Experiment 1: Enzymatic Time Course



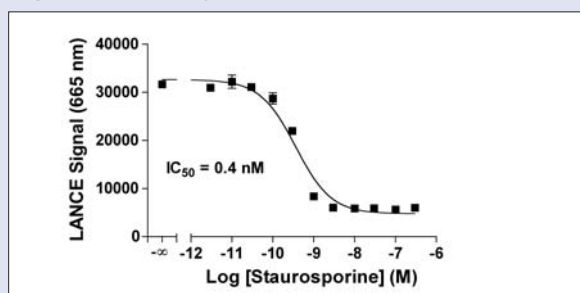
AMPK α 1 enzyme was incubated at concentrations ranging from 0.25 to 4 nM with 50 nM *ULight*-Acetyl-CoA Carboxylase peptide and 20 μ M ATP. Kinase reactions were terminated after 0 to 120 min by the addition of EDTA.

Experiment 2: ATP Titration



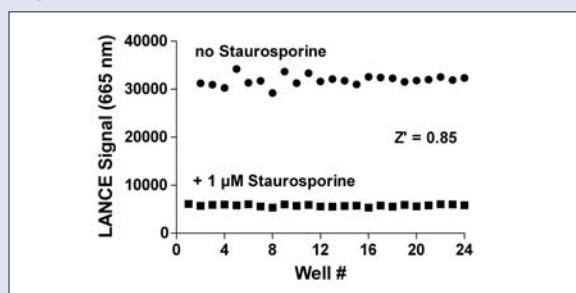
Serial dilutions of ATP ranging from 10 nM to 1 mM were added to 2 nM AMPK α 1 and 50 nM *ULight*-Acetyl-CoA Carboxylase peptide. Kinase reactions were terminated after 30 min by the addition of EDTA.

Experiment 3: Enzyme Inhibition Curve



Serial dilutions of staurosporine ranging from 3 μ M to 300 nM (final concentrations in 2% DMSO) were incubated with 2 nM AMPK α 1, 50 nM *ULight*-Acetyl-CoA Carboxylase peptide and 30 μ M ATP. Kinase reactions were terminated after 30 min by the addition of EDTA.

Experiment 4: Z'-factor Determination



AMPK α 1 enzyme at 2 nM was incubated with 50 nM *ULight*-Acetyl-CoA Carboxylase peptide and 30 μ M ATP with or without 1 μ M staurosporine (final concentrations in 2% DMSO). Kinase reactions were terminated after 30 min by the addition of EDTA.