

U-TRF #11

# LANCE *Ultra* PKC Kinase Assay

LANCE® *Ultra* TR-FRET Technology

PerkinElmer, Inc.  
Montreal, QC  
Canada H3J 1R4

This LANCE *Ultra* kinase assay measures the phosphorylation of a PKC peptide substrate at Ser25.

### Europium-anti-phospho-PKC (Ala25Ser) Peptide Antibody:

- TRF0207-D: 10 µg, 1,562 assay points\*
- TRF0207-M: 100 µg, 15,625 assay points\*

\*40 fmol/assay point

### ULight™-PKC Peptide:

- TRF0108-D: 0.5 nmole, 1,000 assay points\*
- TRF0108-M: 5 nmoles, 10,000 assay points\*

\*0.5 pmol/assay point

### Recognized Motif:

RFARKGpSLRQKNV

Europium-labeled mouse monoclonal antibody recognizing phospho-Ser25 in human PKC in which Ala25 is mutated to Ser and phosphorylated.

### Peptide Sequence:

CRFARKGSLRQKNV

Synthetic peptide derived from amino acids 19-31 of PKC-α; phosphorylation site: Ser25.

### LANCE *Ultra* Kinase Assays:

LANCE *Ultra* time-resolved fluorescence resonance energy transfer (TR-FRET) assays use a proprietary europium chelate donor dye, W1024 (Eu), together with ULight™, a small molecular weight acceptor dye with a red-shifted fluorescent emission.

In this technical note, we present the optimization of a PKC kinase assay using a ULight-labeled peptide substrate. The binding of a Eu-labeled antibody directed against Ser25 phosphorylation of the PKC peptide substrate brings the Eu donor and ULight acceptor dye molecules into close proximity. Upon irradiation 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.

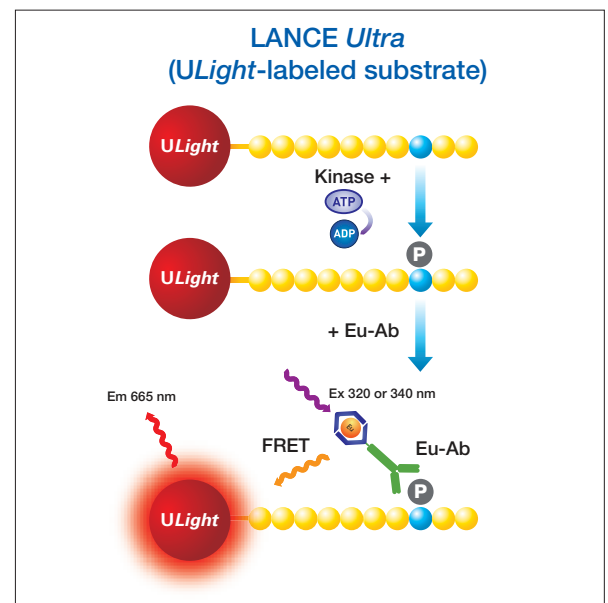


Figure 1. Schematic representation of the LANCE *Ultra* detection of a phosphorylated peptide substrate.

## Development of a PKC Kinase Assay

### Reagents needed for this assay:

Europium-labeled anti-phospho-PKC (Ala25Ser) Peptide Antibody	PerkinElmer # TRF0207
ULight-PKC Peptide	PerkinElmer # TRF0108
PKC- $\alpha$	Invitrogen # P2227
LANCE <sup>®</sup> Detection Buffer, 10X	PerkinElmer # CR97-100
OptiPlate <sup>™</sup> -384, white	PerkinElmer # 6007299
TopSeal <sup>™</sup> -A film	PerkinElmer # 6050195
Kinase Buffer: 50 mM HEPES pH 7.5, 1 mM EGTA, 10 mM MgCl <sub>2</sub> , 2 mM DTT and 0.01% Tween-20.	

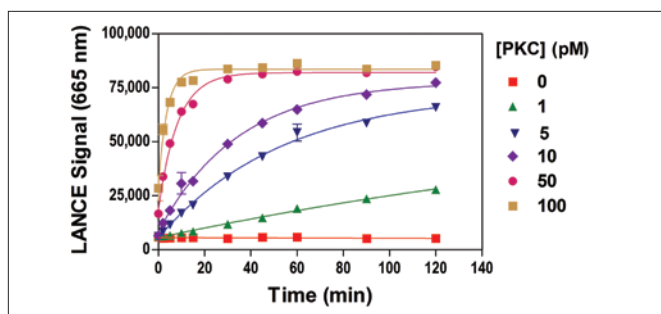
### Standard Protocol

- Dilute the PKC- $\alpha$  kinase, ATP, inhibitors and ULight-PKC peptide in Kinase Buffer.
- Prepare a 4X Detection Mix by diluting the Eu-anti-phospho PKC (Ala25Ser) antibody to 8 nM in 1X LANCE Detection Buffer.

- Add to the wells of a white OptiPlate-384:
  - 5  $\mu$ L of PKC- $\alpha$  enzyme
  - 2.5  $\mu$ L of inhibitor or Kinase Buffer
  - 2.5  $\mu$ L of ULight-PKC substrate/ATP mix (for ATP titration, ATP dilutions are added separately in Kinase Buffer).
- Cover the plate with TopSeal-A film and incubate for 30 min at room temperature (RT).
- Stop kinase reactions by adding 5  $\mu$ L of 40 mM EDTA prepared in 1X LANCE Detection Buffer (Stop Solution). Leave for 5 min at RT.
- Add 5  $\mu$ L of 4X Detection Mix (Eu-anti-phospho-PKC (Ala25Ser) Antibody at a final concentration of 2 nM).
- Cover with TopSeal-A film and incubate for 1 h at RT.
- Remove the TopSeal-A film and read signal with the EnVision<sup>®</sup> Multilabel Reader in TR-FRET mode (excitation at 320 or 340 nm and emission at 665 nm).

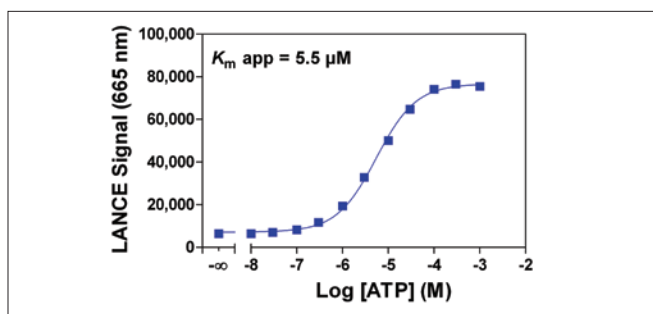
**NOTE:** Eu-labeled antibodies and EDTA can be premixed just before use as a 2X concentrated Stop Solution/Detection Mix to minimize the number of liquid handling steps.

### Experiment 1: Enzymatic Titration and Time Course



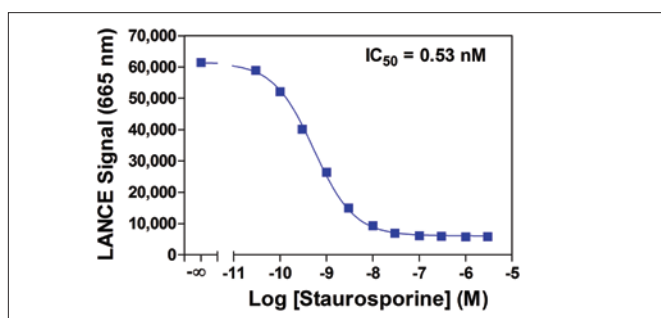
Enzymatic progress curves were produced by incubating PKC- $\alpha$  enzyme at concentrations ranging from 1 to 100 pM with 50 nM ULight-PKC substrate and 10  $\mu$ M ATP. Kinase reactions were terminated at the indicated times by the addition of EDTA. Detection mix was added and signal read after 60 minutes.

### Experiment 2: ATP Titration



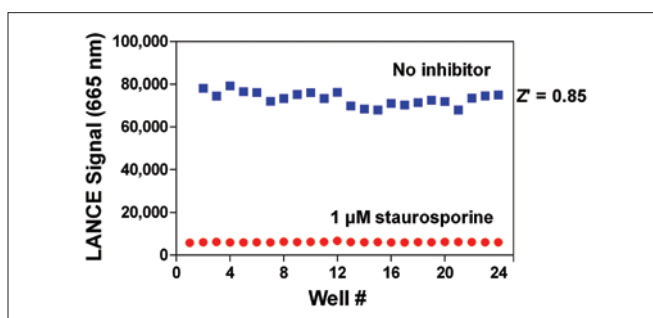
Serial dilutions of ATP ranging from 10 nM to 1 mM were added to 10 pM PKC- $\alpha$  and 50 nM ULight-PKC substrate. Kinase reactions were terminated after 30 min by the addition of EDTA.

### Experiment 3: Enzyme Inhibition Curve



Serial dilutions of staurosporine ranging from 30 pM to 3  $\mu$ M (final concentrations in 1% DMSO) were incubated with 10 pM PKC- $\alpha$ , 50 nM ULight-PKC substrate and 10  $\mu$ M ATP. Kinase reactions were terminated after 30 min by the addition of EDTA.

### Experiment 4: Z'-factor Determination



PKC- $\alpha$  enzyme at 10 pM was incubated with 50 nM ULight-PKC substrate and 10  $\mu$ M ATP, with or without 1  $\mu$ M staurosporine (final concentrations in 1% DMSO). Kinase reactions were terminated after 30 min by the addition of EDTA.