

LANCE *Ultra* Aurora A Kinase Assay

U-TRF #6

LANCE® *Ultra* TR-FRET Technology

PerkinElmer, Inc.

Montreal, QC
Canada H3J 1R4

This LANCE *Ultra* kinase assay measures the phosphorylation of a PLK peptide substrate at Ser137.

Europium-anti-phospho-PLK (Ser137) Antibody:

- TRF0203-D: 10 µg, 1,562 assay points*
- TRF0203-M: 100 µg, 15,562 assay points*
- *40 fmol/assay point
- RECOGNIZED MOTIF: RRRpSLLLE
- Europium-labeled rabbit polyclonal antibody recognizing phospho-Ser137 in human polo-like kinase (PLK)

ULight™-PLK (Ser137) Peptide:

- TRF0110-D: 0.5 nmole, 1,000 assay points*
- TRF0110-M: 5 nmoles, 10,000 assay points*
- *0.5 pmol/assay point
- CORE SEQUENCE MOTIF: RRRSLLE
 - Synthetic peptide containing the residues surrounding Ser137 of polo-like kinase (PLK)
 - Phosphorylation site: Ser137

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 Aurora kinase assay using a ULight-labeled peptide substrate. The binding of an Eu-labeled antibody directed against Ser137 phosphorylation of the PLK 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.

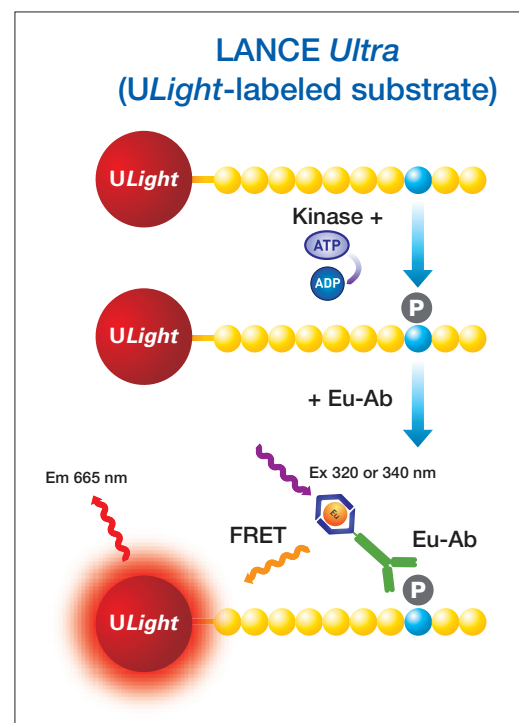


Figure 1. Schematic representation of the LANCE *Ultra* detection of a phosphorylated peptide substrate.

Development of an Aurora A Kinase Assay:

Reagents needed for this assay:

Europium-labeled anti-phospho-PLK (Ser137) Antibody	PerkinElmer # TRF0203
ULight-PLK (Ser137) Peptide	PerkinElmer # TRF0110
Aurora A, active	Millipore # 14-511
LANCE Detection Buffer, 10X	PerkinElmer # CR97-100
OptiPlate™-384, white	PerkinElmer # 6007299
TopSeal-A™ film	PerkinElmer # 6050195

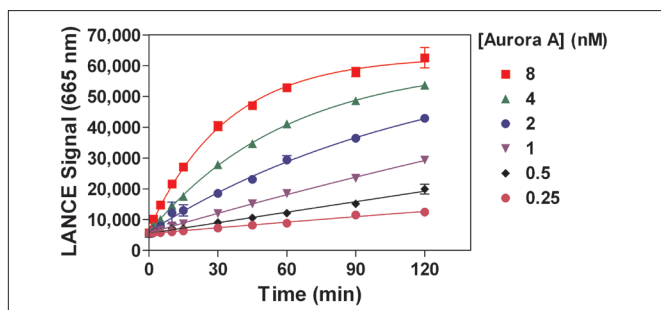
Kinase Buffer: 50 mM HEPES, pH 7.5, 1 mM EGTA, 10 mM MgCl₂, 2 mM DTT and 0.01% Tween-20

Standard Protocol

- Dilute the Aurora A kinase, ATP, inhibitors and ULight-PLK peptide in Kinase Buffer.
- Prepare a 4X Detection Mix by diluting the Eu-anti-phospho-PLK antibody to 8 nM in 1X LANCE Detection Buffer.
- Add to the wells of a white OptiPlate-384:
 - 5 µL of Aurora A enzyme
 - 2.5 µL of inhibitor or Kinase Buffer
 - 2.5 µL of ULight-PLK Peptide/ATP mix (for ATP titration, ATP dilutions are added separately in Kinase Buffer).
- Cover the plate with TopSeal-A film and incubate at room temperature (RT).
- Stop kinase reactions by adding 5 µL of 40 mM EDTA prepared in 1X LANCE Detection Buffer. Leave for 5 min at RT.
- Add 5 µL of Detection Mix (Eu-anti-phospho-PLK antibody at a final concentration of 2 nM).
- Cover with the TopSeal-A film and incubate for 1 h at RT.
- Remove TopSeal-A film and read signal with the EnVision™ Multilabel Reader in TR-FRET mode (excitation at 320 or 340 nm and emission at 665 nm).

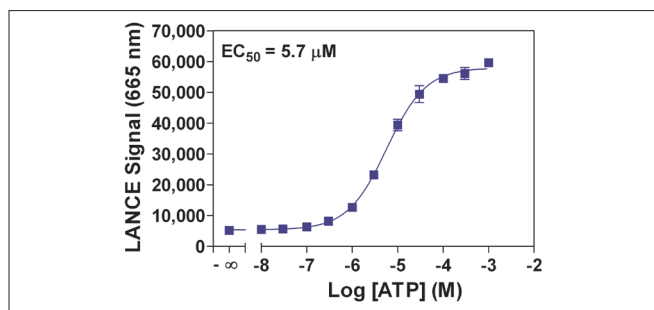
NOTE: Eu-labeled antibodies and EDTA can be premixed just before use as a 2X concentrated Stop/Detection mix to minimize the number of liquid handling steps.

Experiment 1: Enzymatic Titration and Time-Course



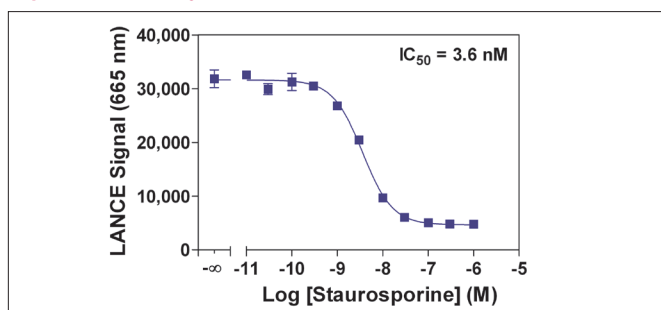
Enzymatic progress curves were produced by incubating Aurora A enzyme at final concentrations ranging from 0.25 to 8 nM with 50 nM ULight-PLK peptide and 20 µM ATP. Kinase reactions were terminated at the indicated times by the addition EDTA. Detection mix was added and signal read after 60 minutes.

Experiment 2: ATP Titration



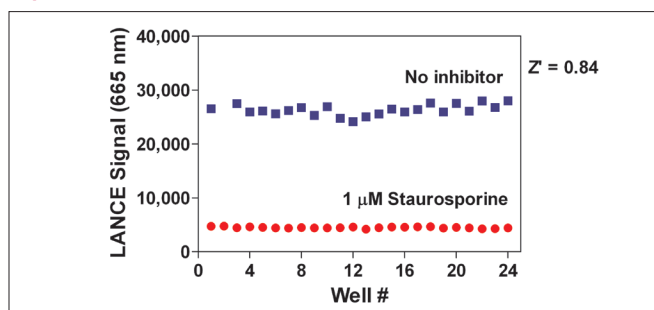
Serial dilutions of ATP ranging from 10 nM to 1 mM were added to 4 nM Aurora A and 50 nM of ULight-PLK peptide. Kinase reactions were terminated after 60 min by the addition of EDTA.

Experiment 3: Enzyme Inhibition



Serial dilutions of staurosporine ranging from 10 pM to 1 µM (final concentrations in 2% DMSO) were incubated with 4 nM Aurora A, 50 nM ULight-PLK peptide and 5 µM ATP. Kinase reactions were terminated after 60 min by the addition of EDTA.

Experiment 4: Z'-factor Determination



Aurora A enzyme at 4 nM was incubated with 50 nM ULight-PLK peptide and 5 µM ATP with or without 1 µM staurosporine (final concentrations in 2% DMSO). Kinase reactions were terminated after 60 min by the addition of EDTA.

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright © 2006-2012, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.