

U-TRF #21

LANCE *Ultra* Haspin Kinase Assay

LANCE® *Ultra* TR-FRET Technology

PerkinElmer, Inc.
Montreal, QC
Canada H3J 1R4

This LANCE *Ultra* kinase assay measures the phosphorylation of a Histone 3 peptide substrate at Thr3.

ULight™-Histone H3 (Thr3/Ser10) Peptide:

- TRF0125-D: 0.5 nmole, 1,000* assay points
- TRF0125-M: 5 nmoles, 10,000* assay points

*0.5 pmol/assay point

Europium-anti-phospho-Histone 3 (Thr3) Antibody:

- TRF0211-D: 10 µg, 1,562* assay points
- TRF0211-M: 100 µg, 15,625* assay points

*40 fmol/assay point

Recognized Motif:

Europium-labeled rabbit monoclonal antibody recognizing phospho-Thr3 in human Histone H3.

Peptide Sequence:

ARTKQTARKSTGGKAPRKQLAGCG

Synthetic peptide containing the residues surrounding Thr3 and Ser10 of human Histone H3; phosphorylation sites: Thr3 and Ser10.

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 Haspin kinase assay using a ULight-labeled peptide substrate. The binding of a Eu-labeled anti-phospho-Histone H3 (Thr3) antibody to the Histone H3 peptide substrate at Thr3 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.

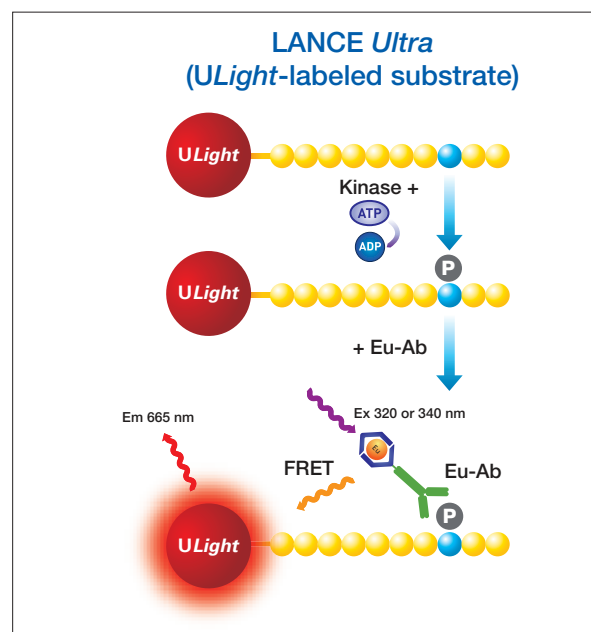


Figure 1. Schematic representation of the LANCE *Ultra* detection of a phosphorylated peptide substrate.

Development of Haspin Kinase Assay

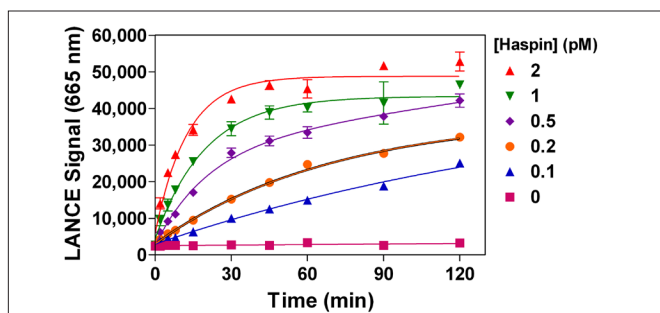
Reagents needed for this assay:

Europium-anti-phospho-Histone H3 (Thr3) Antibody	PerkinElmer # TRF0211
ULight™-Histone H3 (Thr3/Ser10) Peptide	PerkinElmer # TRF0125
Haspin active	Carna # 05-111
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 MgCl ₂ , 2 mM DTT and 0.01% Tween-20	

Standard Protocol

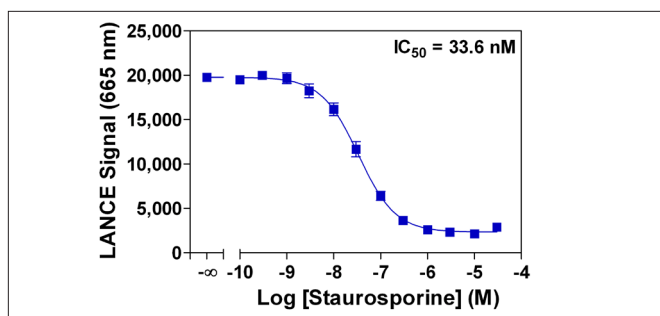
- Dilute the Haspin kinase, ATP, inhibitors and ULight-Histone H3 (Thr3/Ser10) peptide in Kinase Buffer.
- Prepare a 4X Detection Mix by diluting the Eu-anti-phospho-Histone H3 (Thr3) antibody to 8 nM in 1X LANCE Detection Buffer.

Experiment 1: Enzymatic Titration and Time Course



Enzymatic progress curves were produced by incubating Haspin enzyme at concentrations ranging from 0.1 to 2 pM with 50 nM ULight-Histone H3 (Thr3/Ser10) peptide and 100 μM 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 3: Enzyme Inhibition Curve



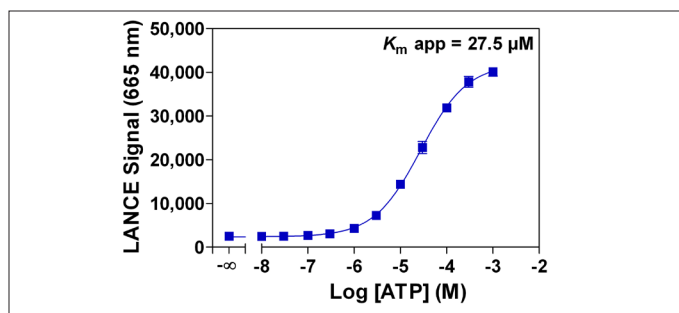
Serial dilutions of staurosporine ranging from 0.1 nM to 30 μM (final concentrations in 1% DMSO) were incubated with 0.5 pM Haspin enzyme, 50 nM ULight-Histone H3 (Thr3/Ser10) peptide and 30 μM ATP. Kinase reactions were terminated after 60 min by the addition of EDTA.

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

- Add to the wells of a white OptiPlate-384:
 - 5 μL of Haspin enzyme
 - 2.5 μL of inhibitor or Kinase Buffer
 - 2.5 μL of ULight-Histone H3 (Thr3/Ser10) 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 24 mM EDTA prepared in 1X LANCE Detection Buffer (Stop Solution). Leave for 5 min at RT.
- Add 5 μL of Detection Mix (Eu-anti-phospho-Histone H3 (Thr3) 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® Multilabel Reader in TR-FRET mode (excitation at 320 or 340 nm and emission at 665 nm).

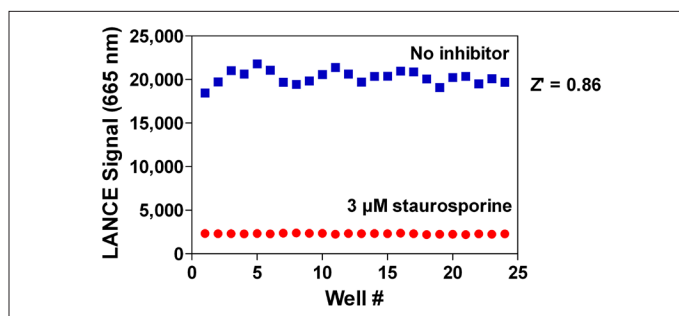
NOTE: Eu-labeled antibodies and EDTA can be premixed just before use as a 2X concentrated Stop Solution/Detection Mix to minimize the number of liquid handling steps.

Experiment 2: ATP Titration



Serial dilutions of ATP ranging from 10 nM to 1 mM were added to 0.5 pM Haspin enzyme and 50 nM ULight-Histone H3 (Thr3/Ser10) peptide. Kinase reactions were terminated after 60 min by the addition of EDTA.

Experiment 4: Z'-factor Determination



Haspin enzyme at 0.5 pM was incubated with 50 nM ULight-Histone H3 (Thr3/Ser10) peptide in kinase assay buffer with 30 μM ATP, and with or without 3 μM staurosporine (final concentrations in 1% DMSO). Kinase reactions were terminated after 60 min by the addition of EDTA.



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

Copyright © 2007-2012, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.