

# LANCE *Ultra* p60<sup>c-src</sup> Kinase Assay

Using *ULight*-Poly GT (4:1) & Europium-Anti-Phospho-Tyrosine (PY20) Antibody

Two LANCE® *Ultra* companion products—two convenient sizes!

*ULight*™-poly GT (4:1):

- **TRF0100-D: 1 nmole, 1,000 assay points\***
- **TRF0100-M: 10 nmoles, 10,000 assay points\***  
\*1 pmol/assay point
- **PEPTIDE SEQUENCE:** [Glu-Tyr (4:1)]<sub>n</sub>
  - Generic substrate for tyrosine kinases
- **VALIDATED FOR KINASES:** p60<sup>c-src</sup>, AXL
- **POTENTIAL SUBSTRATE FOR KINASES:** most tyrosine kinases

Europium-anti-phospho-tyrosine (PY20) antibody:

- **AD0066: 50 µg, 7,800 assay points\***
- **AD0067: 1 mg, 150,000 assay points\***  
\*40 fmol/assay point
- **RECOGNIZED MOTIF: pTyr**
- Mouse monoclonal IgG<sub>2b</sub> antibody directed against phospho-tyrosine.

## 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*, a new innovative small molecular weight acceptor dye with a red-shifted fluorescent emission. In kinase assays, the binding of an Eu-labeled anti-phospho-substrate antibody to phosphorylated *ULight*-labeled substrates brings donor and acceptor molecules into close proximity.

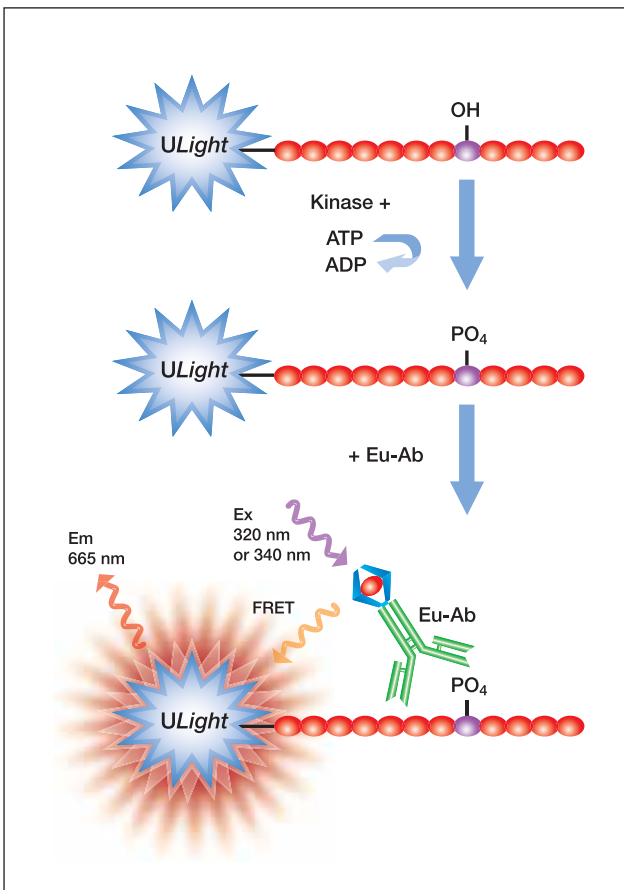
After irradiation of the kinase reaction at 320 nm, the energy from the Eu donor is transferred to the *ULight* acceptor which, in turn, generates light at 665 nm. The intensity of the light emission is proportional to the level of *ULight*-substrate phosphorylation.

## Development of a p60<sup>c-src</sup> Kinase Assay

### Additional Reagents

p60 <sup>c-src</sup> , recombinant	Upstate # 14-326
LANCE Detection Buffer, 10X	PerkinElmer # CR97-100
OptiPlate™-384, white	PerkinElmer # 6007299
TopSeal-A™	PerkinElmer # 6005185

Kinase Buffer: 50 mM Tris-HCl, pH 7.5, 1 mM EGTA, 10 mM MgCl<sub>2</sub>, 2 mM DTT and 0.01% Tween-20



## Suggested Procedure

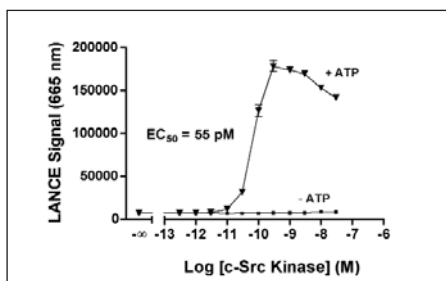
- Dilute kinase, ATP, inhibitors and *ULight*-poly GT (4:1) in Kinase Buffer.
- Dilute antibody (Ab) in LANCE Detection Buffer to 8 nM.
- Add to the wells of a white OptiPlate-384:
  - 5  $\mu$ L of p60<sup>c-src</sup> enzyme,
  - 2.5  $\mu$ L of inhibitor or Kinase Buffer,
  - 2.5  $\mu$ L of *ULight*-poly GT (4:1)/ATP mix (for ATP titration, ATP dilutions are added separately in Kinase Buffer).
- Incubate the enzymatic reaction at room temperature (RT).
- Stop the reaction by adding 5  $\mu$ L of 40 mM EDTA in Detection Buffer. Leave 5 min at RT.
- Add 5  $\mu$ L of the antibody dilution (2 nM final concentration).
- 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).

## Better p60<sup>c-src</sup> Kinase Assays with a Better Technology—LANCE Ultra

For more information about LANCE Ultra, please visit [www.perkinelmer.com/lanceultra](http://www.perkinelmer.com/lanceultra) or contact your local PerkinElmer Sales Representative. Learn more about our comprehensive range of reagents and consumables for drug discovery by visiting [www.perkinelmer.com/drugdiscovery](http://www.perkinelmer.com/drugdiscovery).

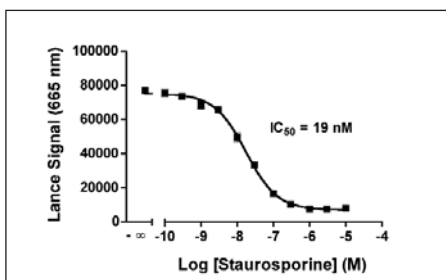
**PerkinElmer Life and Analytical Sciences**  
710 Bridgeport Avenue  
Shelton, CT 06484-4794 USA  
Phone: (800) 762-4000 or  
(+1) 203-925-4602  
[www.perkinelmer.com](http://www.perkinelmer.com)

## Experiment 1: Enzymatic Titration Assay



p60<sup>c-src</sup> enzyme at concentrations ranging from 0.3 pM to 30 nM was incubated with 100 nM *ULight*-poly GT (4:1) and 20  $\mu$ M ATP. Kinase reactions were terminated after 60 min by the addition of EDTA.

## Experiment 3: Enzyme Inhibition Curve

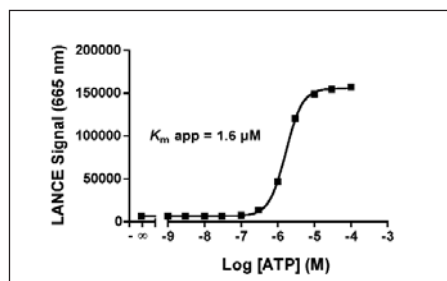


Serial dilutions of staurosporine ranging from 100 pM to 10  $\mu$ M (final concentrations in 2% DMSO) were pre-incubated for 5 min with the p60<sup>c-src</sup> enzyme (300 pM final concentration). Then 100 nM *ULight*-Poly GT (4:1) and 4  $\mu$ M ATP were added. Kinase reactions were terminated after 60 min by the addition of EDTA.

## Complementary Products

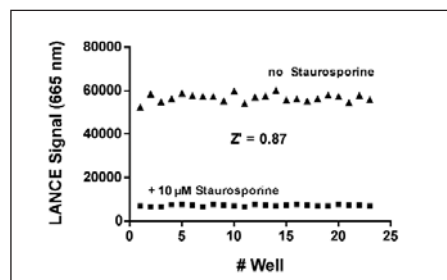
Product	Description	Product No.
Europium-anti-phospho-Tyrosine (PT66)	50 $\mu$ g	AD0068
	1 mg	AD0069
Europium-anti-phospho-Tyrosine (P-Tyr-100)	10 $\mu$ g	AD0203
	50 $\mu$ g	AD0161
	1 mg	AD0162
<i>ULight</i> -poly GAT, [EAY(1:1:1)] <sub>n</sub>	1 nmole	TRF0101-D
	10 nmoles	TRF0101-M

## Experiment 2: ATP Titration



Serial dilutions of ATP ranging from 1 nM to 100  $\mu$ M were added to 300 pM p60<sup>c-src</sup> kinase and 100 nM of *ULight*-poly GT (4:1) substrate. Kinase reactions were terminated after 60 min by the addition of EDTA.

## Experiment 4: Z'-factor Determination:



The p60<sup>c-src</sup> enzyme at 300 pM was incubated with 100 nM *ULight*-poly GT (4:1) substrate in Kinase Assay Buffer with 4  $\mu$ M ATP or 10  $\mu$ M staurosporine and ATP. Reactions were terminated after 60 min by the addition of EDTA.

For a complete listing of our global offices, visit [www.perkinelmer.com/lasoffices](http://www.perkinelmer.com/lasoffices)

©2006 PerkinElmer, Inc. All rights reserved. The PerkinElmer logo and design are registered trademarks of PerkinElmer, Inc. LANCE is a registered trademark and EnVision, OptiPlate, TopSeal-A and *ULight* are trademarks of PerkinElmer, Inc. or its subsidiaries, in the United States and other countries. *ULight* is covered under patent number US2005202565 and US20040166515. All other trademarks not owned by PerkinElmer, Inc. or its subsidiaries that are depicted herein are the property of their respective owners. PerkinElmer reserves the right to change this document at any time without notice and disclaims liability for editorial, pictorial or typographical errors.