

Comparison of LANCE® Ultra with Other TR-FRET Platforms for a Src Tyrosine Kinase Assay

Mireille Legault, Martina Bielefeld-Sévigny, Anne Labonté, Philippe Roby, Ville Laitala, Roger Bossé, Nathalie Rouleau and Lucille Beaudet



1 Introduction

Protein kinases play a major role in normal cellular functions such as cell proliferation, angiogenesis, and cell adhesion. Dysregulation of kinase activity has been shown to be associated with several human diseases including cancer, diabetes and morphological disorders. This makes kinases crucial targets for drug discovery. Analysis of the human genome has revealed the existence of nearly 520 genes encoding kinases. The abundance of these potential therapeutic targets strengthens the need for developing efficient and robust high throughput screening (HTS) platforms for the discovery of kinase modulators.

LANCE® Ultra is PerkinElmer's improved TR-FRET platform enabling the rapid development of robust and sensitive kinase assays. LANCE Ultra reagents include a series of europium (Eu) chelate-labeled anti-phospho-substrate antibodies and several kinase substrates labeled with the new ULight™ acceptor dye. The ULight dye is a small molecular weight fluorescent dye with a red-shifted emission.

This study shows the performance of two LANCE Ultra assay set-ups (ULight-poly GT-based and ULight-streptavidin-based), and compare it to classical LANCE (APC-based), and to an alternative terbium/fluorescein-based TR-FRET platform. Phosphorylation of poly Glu-Tyr (poly GT) by the Src kinase was used as a model system.

2 Materials

Item	Supplier	Cat. Number
➤ ULight™-poly GT	PerkinElmer LAS, Inc.	TRF0100-M
➤ ULight™-streptavidin	PerkinElmer LAS, Inc.	TRF0102-M
➤ SureLight™ APC-SA	PerkinElmer LAS, Inc.	CR130-100
➤ LANCE EU-PY20 Antibody	PerkinElmer LAS, Inc.	AD0066
➤ Biotinylated-poly GT	PerkinElmer LAS, Inc.	custom product
➤ LANCE Detection Buffer 10X	PerkinElmer LAS, Inc.	CR97-100
➤ Fluorescein-Poly GT	Invitrogen Corp.	PV3610
➤ Tb-PY20 antibody	Invitrogen Corp.	PV3552
➤ TR-FRET dilution buffer	Invitrogen Corp.	PV3574
➤ Src, active	Millipore Corp.	14-326
➤ White OptiPlate™-384	PerkinElmer LAS, Inc.	6007290
➤ TopSeal™-A	PerkinElmer LAS, Inc.	6005250
➤ EnVision® MultiLabel Reader	PerkinElmer LAS, Inc.	2103-0010
➤ Mirror: LANCE/DELTA Dual	PerkinElmer LAS, Inc.	2100-4160
➤ Excitation Filter: UV2(TRF) 320 nm	PerkinElmer LAS, Inc.	2100-5060
➤ Emission Filter: Eu 615 nm	PerkinElmer LAS, Inc.	2100-5090
➤ Emission Filter: LANCE 665 nm	PerkinElmer LAS, Inc.	2100-5110

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

3 Methods

LANCE Ultra (ULight-poly GT-based):
In the optimized assay, 1 nM c-Src, 100 nM ULight-poly GT, and ATP (concentrations indicated in text) were pre-diluted in Kinase Buffer and added to a white OptiPlate-384 in a final volume of 10 µL. Plates were incubated for 90 min at RT and kinase reactions were stopped by the addition of 5 µL of EDTA at 40 nM. After 5 min, 5 µL of Eu-labeled PY20 antibody at 8 nM was added for the detection of the phospho-substrate. Detection reactions were incubated for 1 h at RT. The LANCE Ultra signal was detected using the EnVision MultiLabel Reader. Excitation wavelength was set at 320 nm and emission was recorded at 665 nm.

LANCE Ultra (ULight-streptavidin-based):
ULight-streptavidin was used in combination with biotinylated poly GT. Optimized kinase assay conditions were similar to the ones for the assay with ULight-poly GT except that for the capture of the biotinylated substrate. ULight-streptavidin was added at a final concentration of 100 nM together with the Eu-PY20 antibody.

Classical LANCE (APC-streptavidin-based):
SureLight APC-streptavidin was used in combination with biotinylated poly GT. Optimized kinase assay conditions were similar to the ones described above except that for the capture of the biotinylated substrate. SureLight APC-streptavidin was added at a final concentration of 100 nM together with the Eu-PY20 antibody.

TR-FRET (terbium, fluorescein-poly GT-based):
Optimized assay procedures were identical to the ones developed for the LANCE Ultra ULight-poly GT assay except that fluorescein-poly GT and terbium (Tb)-labeled PY20 antibodies were used in place of the ULight substrate and Eu-PY20 antibodies. The fluorescein substrate was diluted in buffer provided by the supplier and Tb-PY20 antibodies were diluted in TR-FRET dilution buffer. The excitation wavelength was set at 320 nm and emission was recorded at 520 nm.

4 LANCE® Ultra Technology

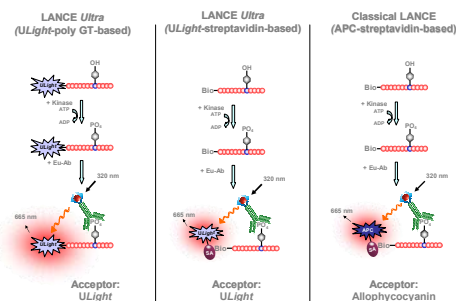
In LANCE TR-FRET assays, the donor dye is an europium-based chelate (Eu chelate). The life-time of the Eu light emission is exceptionally long, allowing for time-delayed measurements. The unique fluorescence properties of Eu chelates make them ideal energy donors in TR-FRET assays.

In classical LANCE assays, the energy emitted upon excitation of an Eu chelate is transferred to an allophycocyanin (APC) acceptor dye molecule which in turn emits light at 665 nm. Although APC allows for the efficient capture and re-emission of the transferred energy, it has some disadvantages. APC is a large protein, whose bulkiness may create steric hindrance in some assay configurations. In addition, small molecules cannot be labeled directly with APC and a bridging assay component, such as streptavidin-APC (APC-SA), must be introduced in the assay set-up.

To overcome these limitations, PerkinElmer has developed the new LANCE Ultra HTS platform where the APC acceptor dye has been replaced by the new ULight dye. This small acceptor dye has spectral characteristics similar to APC and offers the advantage of not being light sensitive. Importantly, its low molecular weight makes it suitable for the direct labeling of molecules of any size. In kinase assays, the ULight dye is coupled either directly to the kinase substrate, or to streptavidin for the capture of a biotinylated substrate. In addition, the red-shifted emission of the ULight dye makes it resistant to color interference which eliminates the need for ratiometric analysis of the data.

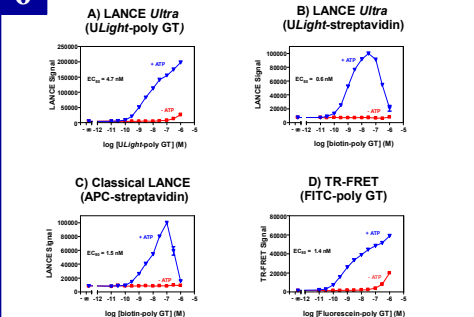
5 Assay Principle

Time-resolved fluorescence resonance energy transfer (TR-FRET) HTS assays are homogeneous proximity assays where the interaction of two dye-labeled binding partners is detected by the energy transfer between a donor and an acceptor dye, and the subsequent light emission by the acceptor dye. Currently, three LANCE TR-FRET platforms are available. They differ principally by the nature of the acceptor dye used for the energy transfer and by the substrate, which can be either directly labeled or biotinylated. Another type of TR-FRET platform currently available uses a terbium chelate as a donor and fluorescein as an acceptor dye.



In LANCE Ultra, ULight-poly GT-based, poly GT is directly labeled with the new ULight acceptor dye (left panel), while in LANCE Ultra, ULight-streptavidin-based, poly GT is biotinylated and captured by a ULight-streptavidin molecule (middle panel). In the classical LANCE APC-based assay, biotinylated poly GT is captured by an APC-SA molecule (right panel). In TR-FRET terbium assays (not shown), peptide substrates are labeled directly with fluorescein. In all four assays, the anti-phosphotyrosine PY20 antibody is labeled with the donor molecule (Eu or Tb chelate). Upon phosphorylation, the poly GT substrate is recognized by the anti-phosphotyrosine antibody bringing the donor and acceptor molecules into close proximity. Excitation of the Eu or Tb chelate leads to an energy transfer to the acceptor molecules resulting in a light emission proportional to the level of substrate phosphorylation.

6 Substrate Titration

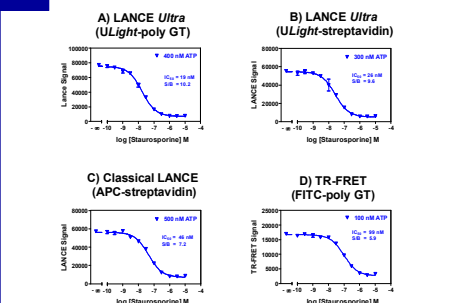


Dilutions of poly GT substrate ranging from 10 pM to 1 µM were incubated in the presence of fixed concentration of c-Src (4 nM) and ATP (20 µM). A) ULight-poly GT, B) and C) biotin-poly GT, D) Fluorescein-poly GT. Detection of the phosphorylated substrates was performed using A) 2 nM of Eu-PY20 antibody, B) 2 nM of Eu-PY20 antibody and 100 nM of ULight-streptavidin, C) 2 nM of Eu-PY20 antibody and 100 nM of APC-streptavidin and D) 2 nM of Tb-PY20 antibody. Note that signal in panels B and C decreases at substrate concentration higher than 100 nM due to the limiting concentration of labeled streptavidin in the assay (hook effect).

EC₅₀ values in the low nanomolar range (0.6 nM to 4.7 nM) were obtained for the poly GT substrate with all four platforms.

Titration of the Src enzyme was also performed for the four assay set-ups (data not shown). Again, similar EC₅₀ values were obtained: 73 pM using ULight-poly GT, 37 pM using ULight-streptavidin, 65 pM using APC-SA and 37 pM using Fluorescein-poly GT.

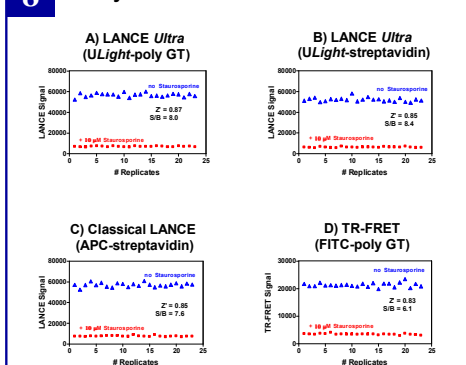
7 Staurosporine Inhibition



Dilutions of staurosporine ranging from 100 pM to 10 µM were pre-incubated for 10 min at RT in the presence of 1 nM of Src kinase. Kinase reactions included 100 nM of poly GT substrate with ATP concentrations at the EC₅₀ value determined for each platform: A) 400 nM, B) 300 nM, C) 500 nM and D) 100 nM. Detection was performed using A) 2 nM of Eu-PY20 antibody, B) 2 nM of Eu-PY20 antibody and 100 nM of ULight-streptavidin, C) 2 nM of Eu-PY20 antibody and 100 nM of APC-streptavidin and D) 2 nM of Tb-PY20 antibody. Signal to background (S/B) ratio is superior in LANCE Ultra assays compared to classical LANCE and TR-FRET assays using a fluorescein-labeled substrate.

IC₅₀ values determined for staurosporine inhibition of the Src kinase were of the same order of magnitude for the four TR-FRET platforms, with values of 19 nM for LANCE Ultra (ULight-poly GT), 26 nM for LANCE Ultra (ULight-streptavidin), 46 nM for classical LANCE, and 99 nM for terbium-based TR-FRET.

8 Assay Performance



Assay performance and robustness were evaluated by calculating the Z' factor¹ with two sets of data: 1) total signal from a kinase reaction incubated in the absence of inhibitor (24 wells) and 2) signal obtained from kinase reactions incubated in the presence of 10 µM staurosporine (24 wells). Src enzyme (1 nM) was pre-incubated in Kinase Buffer in the presence or absence of staurosporine for 10 min at RT. Afterwards, poly GT substrates at a concentration of 100 nM and ATP at concentrations corresponding to the EC₅₀ value (see Figure 7) were then added. Detection was performed using A) 2 nM of Eu-PY20 antibody, B) 2 nM of Eu-PY20 antibody and 100 nM of ULight-streptavidin, C) 2 nM of Eu-PY20 antibody and 100 nM of APC-streptavidin and D) 2 nM of Tb-PY20 antibody.

Z' factors of 0.87 and 0.85 were calculated for the ULight-poly GT assay and ULight-streptavidin assay, respectively. Z' factors obtained using the other two TR-FRET platforms were 0.85 for classical LANCE and 0.83 for fluorescein-poly GT. These results show that excellent assay robustness is achieved with all four TR-FRET platforms.

¹Zhang et al., J Biomol Screen. 1999;4:67-73

9 Conclusions

Optimization of an HTS Src kinase assay using PerkinElmer new LANCE Ultra reagents highlights the robust performance of the new ULight dye as an acceptor for TR-FRET assays.

- The Src assay developed using LANCE Ultra reagents results in performances comparable to classical LANCE, with similar EC₅₀ and IC₅₀ values for poly GT, ATP and staurosporine.
- LANCE Ultra compares favorably to terbium/fluorescein-based TR-FRET platform assays with higher signals, wider assay windows, and an assay robustness demonstrated by Z' factors of 0.85 and above.
- Our data show that, in addition to simplifying assay set-up, the direct labeling of kinase substrates with the smaller ULight dye brings improved assay performance. PerkinElmer also offers ULight custom labeling of proteins, peptides and other small molecules.
- The red-shifted emission properties of the new ULight dye provide minimal screening compound interference. This feature makes LANCE Ultra the platform of choice for HTS kinase assays.
- The new LANCE Ultra reagents offer assay design flexibility with a choice of ULight-labeled kinase substrates and ULight-streptavidin for use with biotinylated substrates.