

# LANCE *Ultra* NEK6 Kinase Assay

## LANCE® *Ultra* TR-FRET Technology

This LANCE *Ultra* kinase assay measures the phosphorylation of a p70 S6K peptide substrate at Thr389.

### **ULight™-p70 S6K (Thr389) Peptide:**

- TRF0126-D: 0.5 nmole, 1,000 assay points\*
- TRF0126-M: 5 nmoles, 10,000 assay points\*

\*0.5 pmol/assay point

### **Europium-anti-phospho-p70 S6K (Thr389) Antibody:**

- TRF0214-D: 10 µg, 1,562 assay points\*
- TRF0214-M: 100 µg, 15,625 assay points\*

\*40 fmol/assay point

### **Recognized Motif:**

FLGF**I**YVAP

Europium-labeled mouse monoclonal antibody recognizing phospho-Thr389 in human p70 S6K.

### **Peptide Motif:**

FLGF**I**YVAP

Synthetic peptide containing the residues surrounding Thr389 of human p70 S6K; phosphorylation site: Thr389.

### **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 NEK6 kinase assay using a *ULight*-labeled peptide substrate. The binding of the Eu-labeled anti-phospho-p70 S6K antibody to the p70 S6K peptide substrate phosphorylated at Thr389 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.

U-TRF #25

PerkinElmer, Inc.  
Montreal, QC  
Canada H3J 1R4

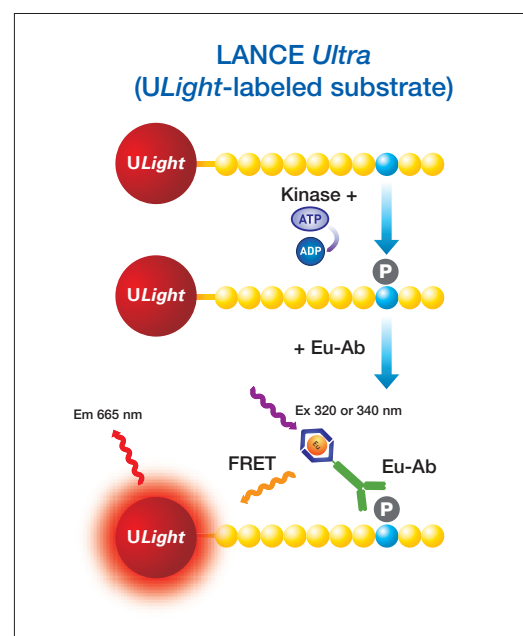


Figure 1. Schematic representation of the LANCE *Ultra* detection of a phosphorylated peptide substrate.

## Development of a NEK6 Kinase Assay

### Reagents needed for this assay:

Europium-labeled anti-phospho-p70 S6K (Thr389) Antibody	PerkinElmer # TRF0214
ULight-p70 S6K (Thr389)	PerkinElmer # TRF0126
NEK6	Carna # 05-130
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 <sub>2</sub> , 2 mM DTT and 0.01% Tween-20.	

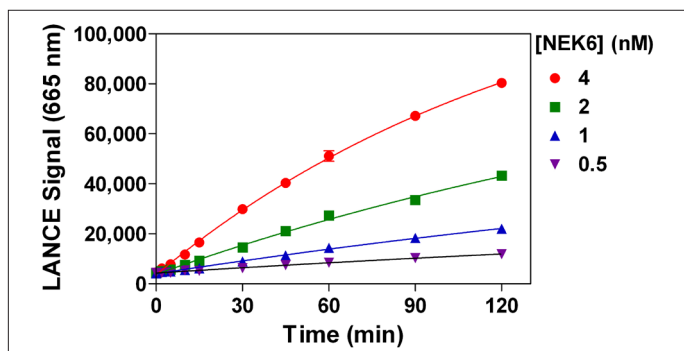
### Standard Protocol

- Dilute the NEK6 enzyme, ATP, inhibitors and ULight-p70 S6K Peptide in Kinase Buffer.
- Prepare a 4X Detection Mix by diluting the Eu-anti-phospho-p70 S6K Antibody to 8 nM in 1X LANCE Detection Buffer.

- Add to the wells of a white OptiPlate-384:
  - 5 µL of NEK6 enzyme
  - 2.5 µL of inhibitor or Kinase Buffer
  - 2.5 µL of ULight-p70 S6K Peptide/ATP mix (for ATP titration, ATP dilutions are added separately in Kinase Buffer).
- Cover the plate with TopSeal-A film and incubate the enzymatic reaction at room temperature (RT).
- Stop kinase reactions by adding 5 µL of 40 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-p70 S6K Antibody at 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 Plate Reader in TR-FRET mode (excitation at 320 or 340 nm and emission at 665 nm).

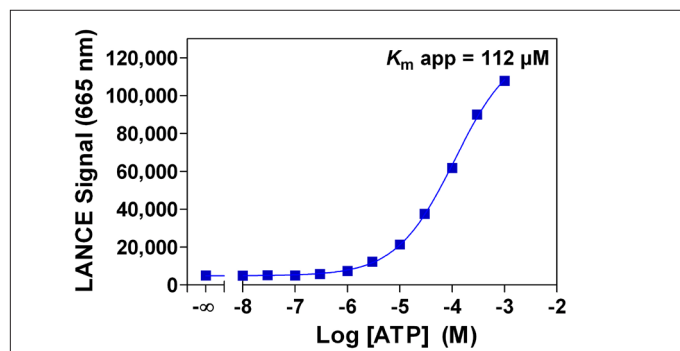
**NOTE:** Eu-labeled antibodies and EDTA can be premixed just before use as a 2X concentrated Stop Solution/Detection Mix to minimize the number of liquid handling steps.

### Experiment 1: Enzymatic Titration and Time Course



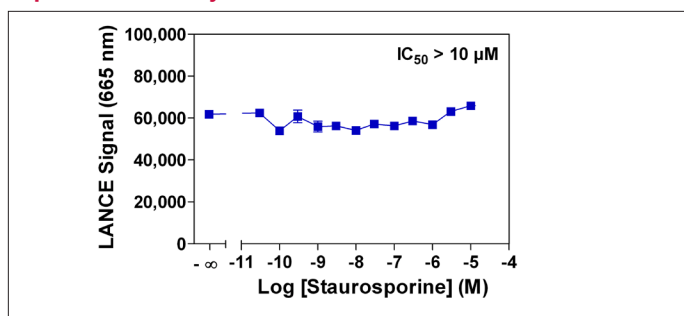
Enzymatic progress curves were produced by incubating NEK6 enzyme at concentrations ranging from 0.5 to 4 nM with 50 nM ULight-p70 S6K Peptide and 100 µM ATP. Kinase reactions were terminated at the indicated times by the addition of EDTA. Detection mix was added and signal read after 60 minutes.

### Experiment 2: ATP Titration



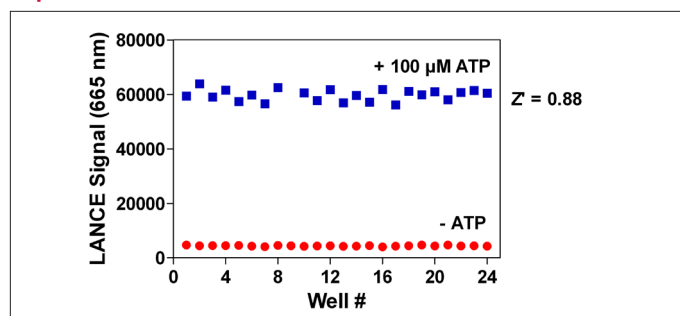
Serial dilutions of ATP ranging from 10 nM to 1 mM were added to 4 nM NEK6 enzyme and 50 nM of ULight-p70 S6K Peptide. Kinase reactions were terminated after 90 min by the addition of EDTA.

### Experiment 3: Enzyme Inhibition Curve



Serial dilutions of staurosporine ranging from 30 pM to 10 µM (final concentrations in 2% DMSO) were incubated with 4 nM NEK6 enzyme, 50 nM ULight-p70 S6K Peptide and 100 µM ATP. Kinase reactions were terminated after 90 min by the addition of EDTA. Staurosporine does not inhibit NEK6 activity at a concentration of 10 µM, consistent with literature data.

### Experiment 4: Z'-factor Determination



NEK6 enzyme at 4 nM was incubated with 50 nM ULight-p70 S6K Peptide in kinase assay buffer with or without 100 µM ATP (final concentrations in 2% DMSO). Kinase reactions were terminated after 90 min by the addition of EDTA.