# **TECHNICAL DATA SHEET**

# THUNDER™ Phospho-EGFR (Y1068) TR-FRET Cell Signaling Assay Kit

CATALOG NUMBERS KIT-EGFRP-100 (100 tests)

KIT-EGFRP-100 (100 tests) KIT-EGFRP-500 (500 tests) KIT-EGFRP-2500 (2500 tests) KIT-EGFRP-5000 (5000 tests) KIT-EGFRP-10000 (10000 tests) Store at -80°C For research use only. Not for use in diagnostic procedures.



# bio**auxilium**

BETTER TOOLS. REAL DISCOVERIES.



#### **PRODUCT DESCRIPTION**

This assay kit measures intracellular levels of **phospho-EGFR (Y1068)** protein in cell lysates using a simple, rapid and sensitive immunoassay based on the homogeneous (no-wash) THUNDER™ TR-FRET technology. The kit is compatible with both adherent and suspension cells.

## SPECIFICITY

This assay kit contains two specific and selective antibodies, one that recognizes **EGFR** phosphorylated at **Tyr1068** and another that recognizes an invariant epitope of **EGFR**.

### SPECIES REACTIVITY

Human (Swiss-Prot Acc. P00533 and Entrez-Gene Id 1956).

Other species should be tested on a case-by-case basis.

## **TR-FRET ASSAY PRINCIPLE**

The Phospho-EGFR (Y1068) assay kit is a homogeneous timeresolved Förster resonance energy transfer (TR-FRET) sandwich immunoassay (Figure 1). The THUNDER™ Cell Signaling assay workflow consists of 3 steps (Figure 2). Following cell treatment, cells are first lysed with the specific Lysis Buffer provided in the kit. Then Phospho-EGFR (Y1068) in the cell lysates is detected with a pair of fluorophore-labeled antibodies in a simple "add-incubatemeasure" format (single-step reagent addition; no wash steps). One antibody is labeled with a donor fluorophore (Europium chelate: Eu-Abl) and the second with a far-red acceptor fluorophore (FR-Ab2). The binding of the two labeled antibodies to distinct epitopes on the target protein takes place in solution and brings the two dyes into close proximity. Excitation of the donor Europium chelate molecules with a flash lamp (320 or 340 nm) or a laser (337 nm) triggers a FRET from the donor to the acceptor molecules, which in turn emit a TR-FRET signal at 665 nm. Residual energy from the Eu chelate generates light at 615 nm. The signal at 665 nm is proportional to the concentration of Phospho-EGFR (Y1068) in the cell lysate. Data can be expressed as either the signal at 665 nm or the 665 nm/615 nm ratio.

STEP 1	STEP 2 STEP 3	
Cell treatment	Cell lysis	Protein detection
<ul> <li>Seed adherent cells in culture plate</li> <li>Add media +/- compound</li> <li>Incubate for optimized time</li> </ul>	Remove media     Add 1X     Supplemented     Lysis Buffer 1     Incubate for 30 min	<ul> <li>Transfer lysate         <ul> <li>(15 μL) to detection plate</li> <li>Add 4X Antibody Mix (5 μL)</li> <li>Incubate for 1 h</li> <li>Read TR-FRET signal</li> </ul> </li> </ul>

Figure 2 Assay workflow using the 2-plate (transfer) protocol.

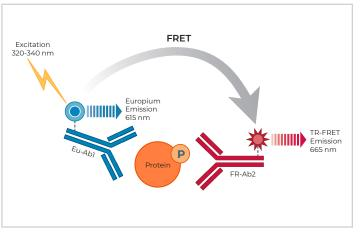


Figure 1 Schematic representation of the TR-FRET cell signaling assay principle.

KIT COMPONENTS	100 points*	500 points*	
Eu-labeled Phospho-EGFR (Y1068) antibody (Eu-Ab1)	5 µL	25 µL	
Acceptor-labeled Phospho-EGFR (Y1068) antibody (FR-Ab2)	20 µL	100 µL	
Lysis Buffer 1 (5X)	lmL	5 mL	
Detection Buffer (10X)	50 µL	250 µL	
Positive control cell lysate	100 µL	200 µL	
Phosphatase Inhibitor Cocktail (100X)	50 µL	250 µL	

\* The number of assay points is based on an assay volume of 20  $\mu$ L in half-area 96-well or low-volume 384-well assay plates using the kit components at the recommended concentrations (refer to the User Manual).

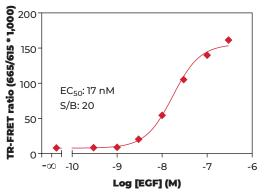
# **TECHNICAL DATA SHEET** Phospho-EGFR (Y1068)

# **VALIDATION DATA**

This assay kit has been validated for the relative quantification of phospho-EGFR (Y1068) in A431 cell lysates using the 2-plate assay protocol.

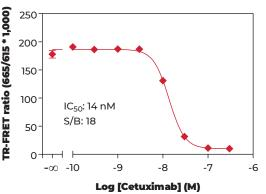
- · Adherent cells were cultured overnight in a 96-well tissue culture plate (DMEM +10% FBS).
- · Following cell treatment, the media was removed and cells were lysed with the **1X Lysis Buffer 1** (50  $\mu$ L) supplemented with the 100X Phosphatase Inhibitor Cocktail diluted at 1X.
- · Following a 30-min incubation at room temperature (RT) on an orbital shaker (400 rpm), lysates (15  $\mu$ L) were then transferred to a 384-well assay plate followed by addition of the labeled antibodies Eu-Abl and FR-Ab2 (5 µL) for detection of phospho-EGFR (Y1068).
- The plate was incubated at RT for **1 hour** and the TR-FRET signal was recorded at 665 and 615 nm (EnVision®; lamp excitation).

# STIMULATION OF PHOSPHO-EGFR (Y1068) IN A431 CELLS

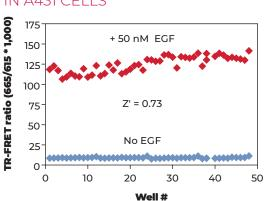


A431 cells (40,000 cells/well; in triplicate) were incubated with serial dilutions of EGF for 10 min at RT. Data show that treatment of A431 cells with EGF stimulates phosphorylation of EGFR at Y1068.

# INHIBITION OF PHOSPHO-EGFR (Y1068) IN A431 CELLS



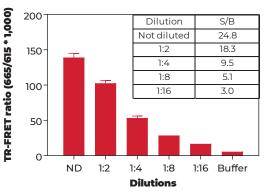
A431 cells (40,000 cells/well; in triplicate) were incubated with serial dilutions of Cetuximab for 30 min at RT. Cells were then stimulated with 50 nM of EGF for 10 min at RT. Data show that treatment of A431 cells with Cetuximab inhibits phosphorylation of EGFR at Y1068 by EGF.



Z'-FACTOR DETERMINATION IN A431 CELLS

A431 cells (40,000 cells/well; in triplicate) were incubated without or with 50 nM of EGF for 10 min at RT. The Z' factor value was determined using a total of 48 wells for each treatment group. The Z'-factor value of 0.73 indicates that the assay is robust and suitable for HTS.

# A431 CONTROL LYSATE TITRATION (OC TEST)



Quality Control: the Phospho-EGFR (Y1068) assay kit is routinely tested against EGF-treated A431 lysates. A431 cells were cultured in a T175 flask to 100% confluence and stimulated with 100 nM of EGF for 10 min at RT. Following cell lysis using 4 mL of 1X Lysis Buffer 1, lysates were serially diluted with 1X Lysis Buffer 1 and tested in triplicate. Data show a linear relationship between lysate dilutions and TR-FRET ratio values.



#### FOR MORE INFORMATION ON DEVELOPING AND OPTIMIZING TR-FRET CELL SIGNALING ASSAYS. CONSULT THE USER MANUAL.

7171 Frederick-Banting, Suite 2115, Montréal (Québec) H4S 1Z9 CANADA bioauxilium.com 1-866-610-6933 x 103