

## TECHNICAL DATA SHEET



bioauxilium

BETTER TOOLS. REAL DISCOVERIES.

# THUNDER™ Phospho-STAT5 (Y694/Y699) + Total STAT5 TR-FRET Cell Signaling Assay Kit

**CATALOG NUMBERS** KIT-STAT5PT-500  
400 points for phospho-STAT5 and 100 points for total STAT5

Store at -80°C  
For research use only.  
Not for use in diagnostic procedures.

## PRODUCT DESCRIPTION

This assay kit measures intracellular levels of **Phospho-STAT5 (Y694/Y699)** and **total STAT5** 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 antibody pairs, one that recognizes **STAT5** phosphorylated at **Tyr694** and **Tyr699** and another that recognizes **total** (both phosphorylated and unphosphorylated) **STAT5**.

## SPECIES REACTIVITY

Human, Mouse (Swiss-Prot Acc.: P42229, P51692; Entrez-Gene Id: 6776, 6777).

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

## TR-FRET ASSAY PRINCIPLE

The **Phospho-STAT5 (Y694/Y699) + Total STAT5** assay kit is a homogeneous time-resolved 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-STAT5 (Y694/Y699)** and **Total STAT5** in the cell lysates are detected in separate wells with two pairs of fluorophore-labeled antibodies in a simple "add-incubate-measure" format (single-step reagent addition; no wash steps). For detection of the phosphorylated protein, one antibody is labeled with a donor fluorophore (Europium chelate; Eu-Ab1) and the second with a far-red acceptor fluorophore (FR-Ab2). The same approach is used for the second antibody pair detecting the total protein (Eu-Ab3 and FR-Ab4). The binding of the two matched labeled antibodies to distinct epitopes on the target protein (either **phospho-STAT5** or **total STAT5**) 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-STAT5 (Y694/Y699)** and **Total STAT5** in the cell lysate. Data can be expressed as either the signal at 665 nm or the 665 nm/615 nm ratio.

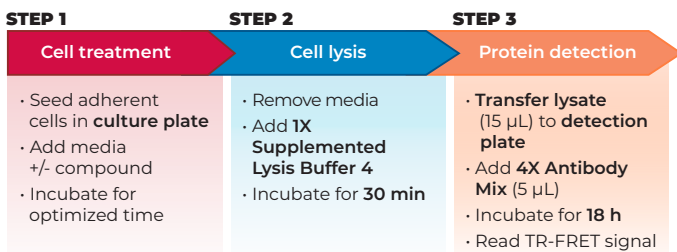


Figure 2 Assay workflow using the 2-plate (transfer) protocol.

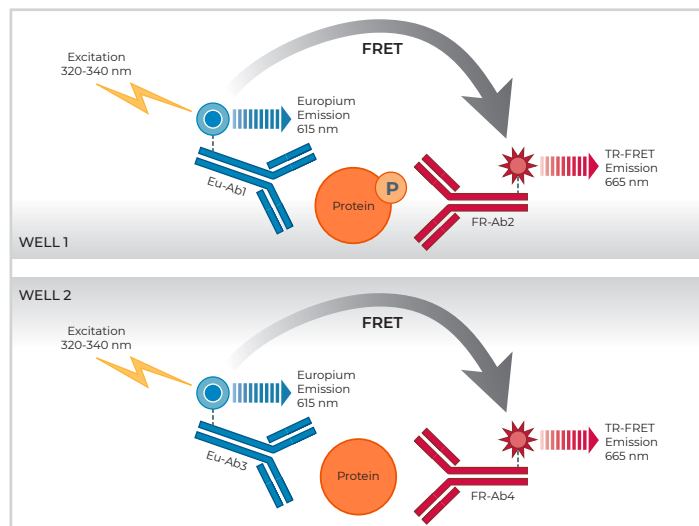


Fig. 1 Schematic representation of the TR-FRET cell signaling assay principle.

## KIT COMPONENTS

	500 points*
Eu-labeled Phospho-STAT5 (Y694/Y699) antibody (Eu-Ab1)	20 µL
Acceptor-labeled Phospho-STAT5 (Y694/Y699) antibody (FR-Ab2)	80 µL
Eu-labeled total-STAT5 antibody (Eu-Ab3)	5 µL
Acceptor-labeled total-STAT5 antibody (FR-Ab4)	20 µL
Lysis Buffer 4 (5X)	5 mL
Detection Buffer (10X)	250 µL
Positive control cell lysate	200 µL
Phosphatase Inhibitor Cocktail (100X)	250 µL

\* The number of assay points is based on an assay volume of 20 µL 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-STAT5 (Y694/Y699) + Total STAT5

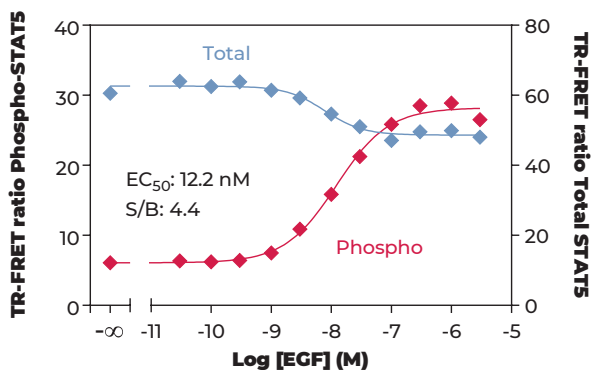
## VALIDATION DATA

This assay kit has been validated for the relative quantification of phospho-STAT5 (Y694/Y699) and total STAT5 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 4** (25  $\mu$ L) supplemented with the phosphatase inhibitors sodium fluoride (1 mM) and sodium orthovanadate (2 mM).
- Following a **30-min** incubation at room temperature (RT) on an

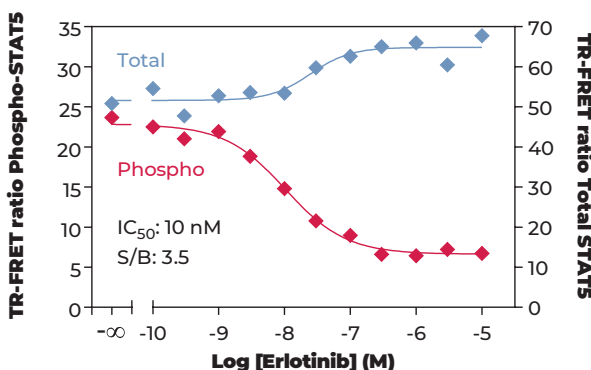
- orbital shaker (400 rpm), lysates (15  $\mu$ L) were then transferred to a 384-well assay plate followed by addition to separate wells of either the labeled antibodies Eu-Ab1 and FR-Ab2 (5  $\mu$ L) for detection of phospho-STAT5 (Y694/Y699) or Eu-Ab3 and FR-Ab4 (5  $\mu$ L) for detection of total STAT5.
- The plate was incubated at RT for **18 hours** and the TR-FRET signal was recorded at 665 and 615 nm (EnVision<sup>®</sup>; lamp excitation).

## STIMULATION OF PHOSPHO-STAT5 (Y694/Y699) IN A431 CELLS



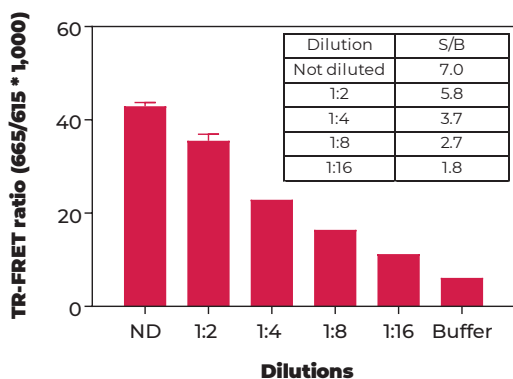
A431 cells (75,000 cells/well; in triplicate) were incubated with serial dilutions of EGF for 10 min at RT. Data show that treatment of A431 cells with EGF stimulates phosphorylation of STAT5 at Y694/Y699, but does not have a major effect on the levels of total STAT5.

## INHIBITION OF PHOSPHO-STAT5 (Y694/Y699) IN A431 CELLS



A431 cells (75,000 cells/well; in triplicate) were incubated with serial dilutions of the inhibitor Erlotinib for 15 min at RT. Cells were then stimulated with 73 nM of EGF for 10 min at RT. Data show that treatment of A431 cells with Erlotinib inhibits phosphorylation of STAT5 at Y694/Y699 by EGF, but does not have a major effect on the levels of total STAT5.

## A431 CONTROL LYSATE TITRATION (QC TEST) PHOSPHO-STAT5 (Y694/Y699)



Quality Control: the Phospho-STAT5 (Y694/Y699) + Total STAT5 assay kit is routinely tested against EGF-treated A431 lysates. A431 cells were cultured in a T175 flask to 100% confluence and stimulated with 300 nM of EGF for 10 min at RT. Following cell lysis using 4 mL of 1X Lysis Buffer 4, lysates were serially diluted with 1X Lysis Buffer 4 and tested in triplicate and in separate wells for phospho-STAT5 (Y694/Y699) and total STAT5. Data show a linear relationship between lysate dilutions and TR-FRET ratio values. Note that due to the very high sensitivity of the Total STAT5 kit, lysates from the T175 flask required at least a 1:2 pre-dilution in order to be within the dynamic assay range.

## A431 CONTROL LYSATE TITRATION (QC TEST) TOTAL STAT5

