

## TECHNICAL DATA SHEET



bioauxilium

BETTER TOOLS. REAL DISCOVERIES.

# THUNDER™ Total STAT3 TR-FRET Cell Signaling Assay Kit

**CATALOG NUMBERS** KIT-STAT3T-100 (100 tests)  
KIT-STAT3T-500 (500 tests)

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

## PRODUCT DESCRIPTION

This assay kit measures intracellular levels of **total STAT3** 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 that recognize **total** (both phosphorylated and unphosphorylated) **STAT3**.

## SPECIES REACTIVITY

Human (Swiss-Prot Acc.: P40763; Entrez-Gene Id: 6774).

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

## TR-FRET ASSAY PRINCIPLE

The **Total STAT3** 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 **Total STAT3** in the cell lysates is detected with a pair of fluorophore-labeled antibodies in a simple "add-incubate-measure" format (single-step reagent addition; no wash steps). One antibody is labeled with a donor fluorophore (Europium chelate; Eu-Ab1) 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 **Total STAT3** in the cell lysate. Data can be expressed as either the signal at 665 nm or the 665 nm/615 nm ratio.

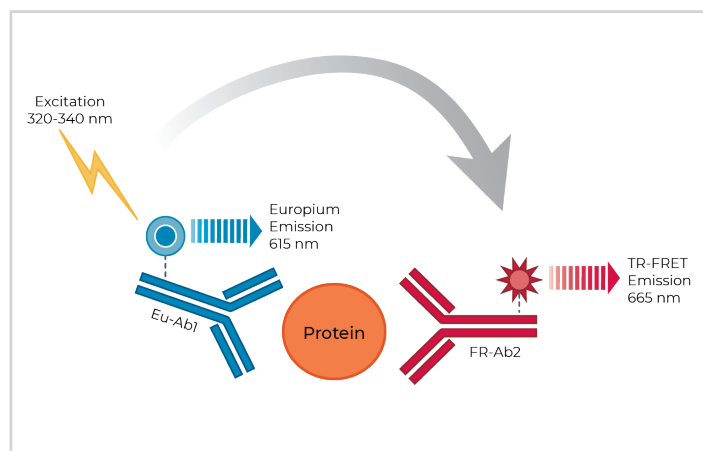


Figure 1 Schematic representation of the TR-FRET cell signaling assay principle.

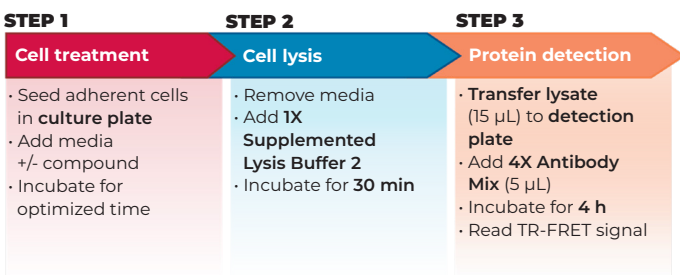


Figure 2 Assay workflow using the 2-plate (transfer) protocol.

## KIT COMPONENTS

	100 points*	500 points*
Eu-labeled total-STAT3 antibody (Eu-Ab1)	5 µL	25 µL
Acceptor-labeled total-STAT3 antibody (FR-Ab2)	20 µL	100 µL
Lysis Buffer 2 (5X)	1 mL	5 mL
Detection Buffer (10X)	50 µL	250 µL
Positive control cell lysate	100 µL	500 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).

## VALIDATION DATA

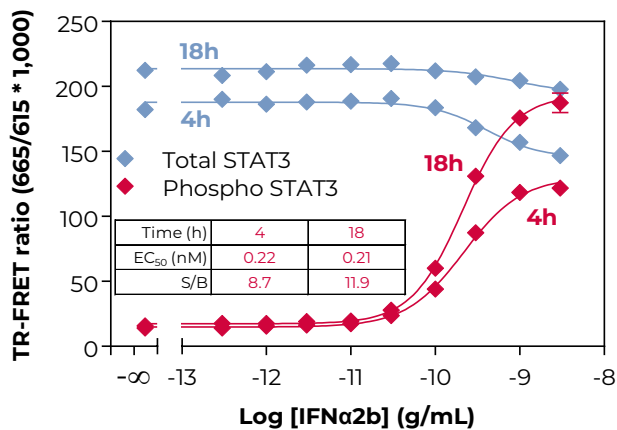
This assay kit has been validated for the relative quantification of total STAT3 in HeLa cell lysates using the 2 plate assay protocol.

- 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 2** (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-Ab1 and FR-Ab2 (5  $\mu$ L) for detection of total STAT3.

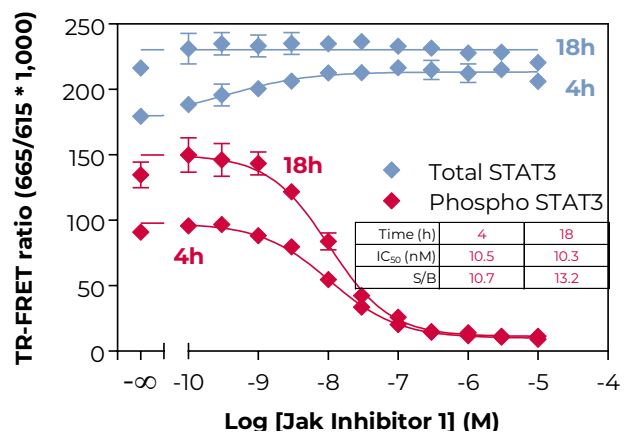
- The plate was incubated at RT for **4 and 18 hours** and the TR-FRET signal was recorded at 665 and 615 nm (EnVision<sup>®</sup>; laser excitation).

## STIMULATION OF PHOSPHO-STAT3 (Y705) IN HELA CELLS



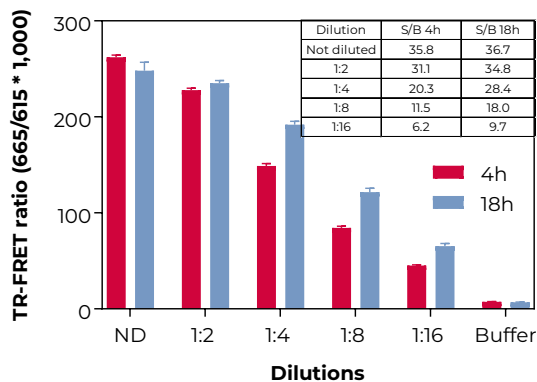
HeLa cells (80,000 cells/well; in triplicate) were incubated with serial dilutions of IFN $\alpha$  for 60 min at RT. Data show that treatment of HeLa cells with IFN $\alpha$  stimulates phosphorylation of STAT3 (Y705), but does not affect the levels of total STAT3.

## INHIBITION OF PHOSPHO-STAT3 (Y705) IN HELA CELLS



HeLa cells (80,000 cells/well; in triplicate) were incubated with serial dilutions of JAK Inhibitor 1 for 30 min at RT. Cells were then stimulated with 0.8 nM of IFN $\alpha$  for 60 min at RT. Data show that treatment of HeLa cells with JAK Inhibitor 1 inhibits phosphorylation of STAT3 (Y705) by IFN $\alpha$ , but does not affect the levels of total STAT3.

## HELA CONTROL LYSATE TITRATION (QC TEST)



Quality Control: the Total STAT3 assay kit is routinely tested against IFN $\alpha$ -treated HeLa lysates. HeLa cells were cultured in a T175 flask to 90% confluence and stimulated with 3 nM of IFN $\alpha$  for 60 min at RT. Following cell lysis using 4 mL of 1X Lysis Buffer 2, lysates were serially diluted with 1X Lysis Buffer 2 and tested in triplicate and in separate wells for phospho-STAT3 (Y705) and total STAT3. Data show a linear relationship between lysate dilutions and TR-FRET ratio values. Note that due to the very high sensitivity of the Total STAT3 kit, lysates from the T175 flask required at least a 1:4 pre-dilution in order to be within the dynamic assay range.

