

## TECHNICAL DATA SHEET

# THUNDER™ Total 4EBP1 TR-FRET Cell Signaling Assay Kit



bioauxilium  
BETTER TOOLS. REAL DISCOVERIES.

**CATALOG NUMBERS** KIT-4EBP1T-100 (100 tests)  
KIT-4EBP1T-500 (500 tests)  
KIT-4EBP1T-2500 (2500 tests)  
KIT-4EBP1T-5000 (5000 tests)  
KIT-4EBP1T-10000 (10000 tests)

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

[Visit product page](#)

## PRODUCT DESCRIPTION

This assay kit measures intracellular levels of **total 4EBP1** 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

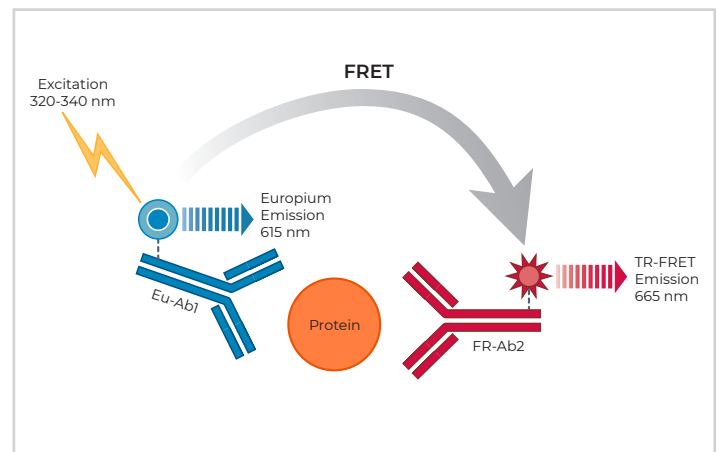
This assay kit contains two specific and selective antibodies that recognize **total** (both phosphorylated and unphosphorylated) **4EBP1**.

## SPECIES REACTIVITY

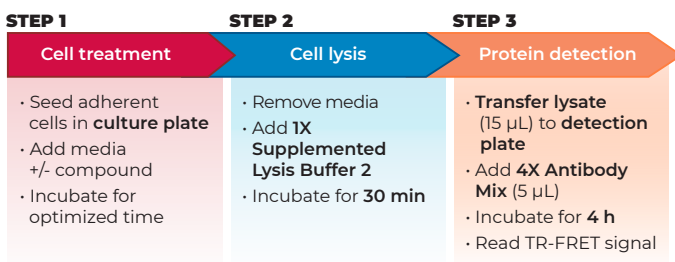
Human (Swiss-Prot Acc.: Q13541; Entrez-Gene Id: 1978).  
Other species should be tested on a case-by-case basis.

## TR-FRET ASSAY PRINCIPLE

The **Total 4EBP1** 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 4EBP1** 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 4EBP1** 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-4EBP1 antibody (Eu-Ab1)       | 5 µL        | 25 µL       |
| Acceptor-labeled total-4EBP1 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      | 200 µL      |
| Phosphatase Inhibitor Cocktail (100X)          | 50 µL       | 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).

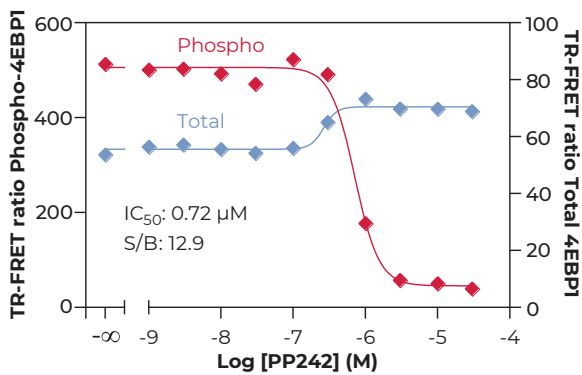
Visit product page

VALIDATION DATA

This assay kit has been validated for the relative quantification of total 4EBP1 in HEK293 and A431 cell lysates using the 2 plate assay protocol.

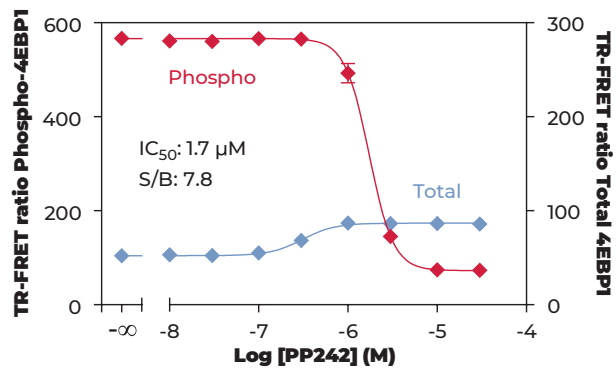
- Adherent cells were cultured overnight in a 96-well tissue culture plate (EMEM +10% FBS for HEK293; DMEM +10% FBS for A431). The plates used for HEK293 were coated with poly-L-lysine.
- Following cell treatment, the media was removed and cells were lysed with the 1X **Lysis Buffer 2** (50  $\mu$ L) 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 of the labeled antibodies Eu-Ab1 and FR-Ab2 (5  $\mu$ L) for detection of total 4EBP1.
- The plate was incubated at RT for **4 hours** and the TR-FRET signal was recorded at 665 and 615 nm (EnVision<sup>®</sup>; lamp excitation).

INHIBITION OF PHOSPHO-4EBP1 (T37/T46) IN A431 CELLS



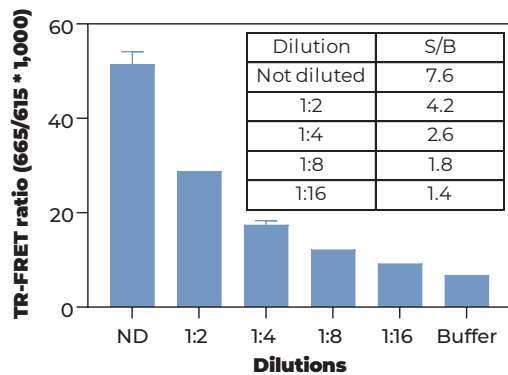
A431 cells (25,000 cells/well; in triplicate) were incubated with serial dilutions of PP242 for 3 hours at 37°C. Data show that treatment of A431 cells with PP242 inhibits phosphorylation of 4EBP1 at T37/T46, but does not have a major effect on the levels of total 4EBP1.

INHIBITION OF PHOSPHO-4EBP1 (T37/T46) IN HEK293 CELLS



HEK293 cells (60,000 cells/well; in triplicate) were incubated with serial dilutions of PP242 for 3 hours at 37°C. Data show that treatment of HEK293 cells with PP242 inhibits phosphorylation of 4EBP1 at T37/T46, but does not have a major effect on the levels of total 4EBP1.

A431 CONTROL LYSATE TITRATION (QC TEST)



Quality Control: the Total 4EBP1 assay kit is routinely tested against A431 lysates. A431 cells were cultured in a T175 flask to 70% confluence. Following cell lysis using 4 mL of 1X Lysis Buffer 2, lysates were serially diluted with 1X Lysis Buffer 2 and tested in triplicate. Data show a linear relationship between lysate dilutions and TR-FRET ratio values.

