

## TECHNICAL DATA SHEET

# THUNDER™ Phospho-JNK1/2/3 (T183/Y185) + Total JNK1/2/3 TR-FRET Cell Signaling Assay Kit



bioauxilium  
BETTER TOOLS. REAL DISCOVERIES.

**CATALOG NUMBERS** KIT-JNKPT-500  
400 points for phospho-JNK and 100 points for total JNK

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

## PRODUCT DESCRIPTION

This assay kit measures intracellular levels of **phospho-JNK1/2/3 (T183/Y185)** and **total JNK1/2/3** 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 **JNK1/2/3** phosphorylated at **Thr183** and **Tyr185** and another that recognizes **total** (both phosphorylated and unphosphorylated) **JNK1/2/3**.

## SPECIES REACTIVITY

Human (Swiss-Prot Acc.: P45983, P45984, P53779; Entrez-Gene Id: 5599, 5601, 5602).

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

## TR-FRET ASSAY PRINCIPLE

The **Phospho-JNK1/2/3 (T183/Y185) + Total JNK1/2/3** 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-JNK1/2/3 (T183/Y185)** and **Total JNK1/2/3** 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-JNK1/2/3** or **total JNK1/2/3**) 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-JNK1/2/3 (T183/Y185)** and **Total JNK1/2/3** in the cell lysate. Data can be expressed as either the signal at 665 nm or the 665 nm/615 nm ratio.

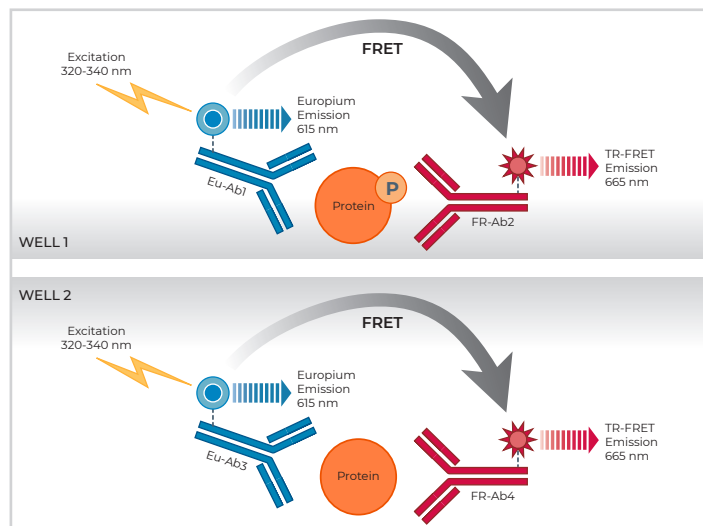


Fig. 1 Schematic representation of the TR-FRET cell signaling assay principle.

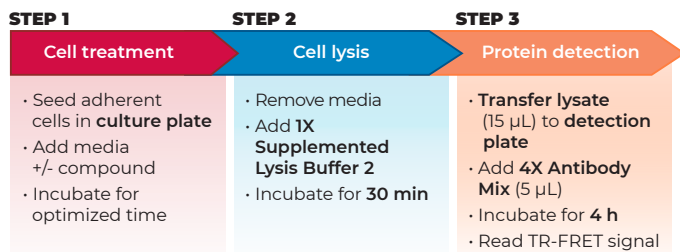


Figure 2 Assay workflow using the 2-plate (transfer) protocol.

## KIT COMPONENTS

	500 points*
Eu-labeled phospho-JNK1/2/3 (T183/Y185) antibody (Eu-Ab1)	20 µL
Acceptor-labeled phospho-JNK1/2/3 (T183/Y185) antibody (FR-Ab2)	80 µL
Eu-labeled total-JNK1/2/3 antibody (Eu-Ab3)	5 µL
Acceptor-labeled total-JNK1/2/3 antibody (FR-Ab4)	20 µL
Lysis Buffer 2 (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-JNK1/2/3 (T183/Y185) + Total JNK1/2/3

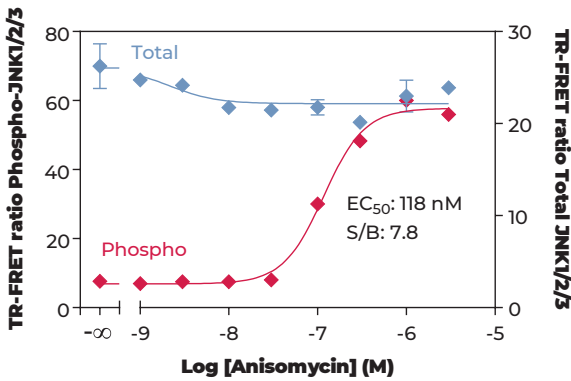
## VALIDATION DATA

This assay kit has been validated for the relative quantification of phospho-JNK1/2/3 (T183/Y185) and total JNK1/2/3 in HEK293 cell lysates using the 2-plate assay protocol.

- Adherent cells were cultured overnight in a 96-well tissue culture plate coated with poly-L-lysine (EMEM +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 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

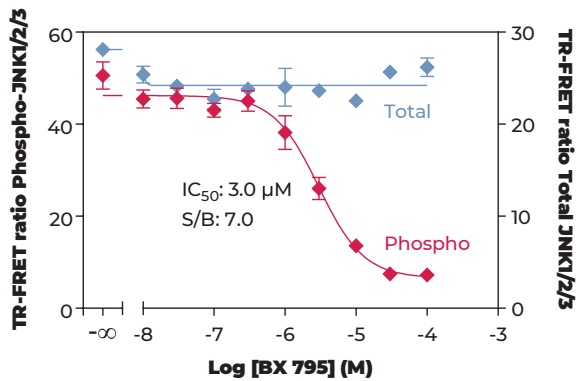
- to a 384-well assay plate followed by addition to separate wells of either the labeled antibodies Eu-Ab1 and FR-Ab2 (5  $\mu$ L) for detection of phospho-JNK1/2/3 (T183/Y185) or Eu-Ab3 and FR-Ab4 (5  $\mu$ L) for detection of total JNK1/2/3.
- The plate was incubated at RT for **4 hours** and the TR-FRET signal was recorded at 665 and 615 nm (EnVision<sup>®</sup>; lamp excitation).

## STIMULATION OF PHOSPHO-JNK1/2/3 (T183/Y185) IN HEK293 CELLS



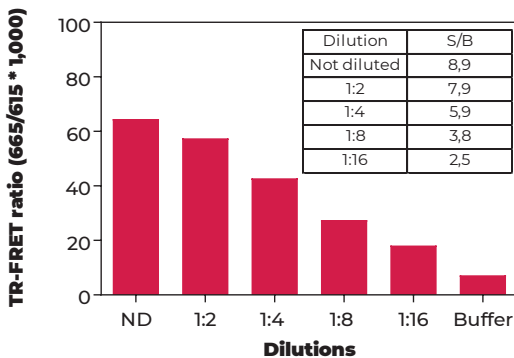
HEK293 cells (50,000 cells/well; in triplicate) were incubated with serial dilutions of Anisomycin for 60 min at RT. Data show that treatment of HEK293 cells with Anisomycin stimulates phosphorylation of JNK1/2/3 at T183/Y185 but does not affect the levels of total JNK1/2/3.

## INHIBITION OF PHOSPHO-JNK1/2/3 (T183/Y185) IN HEK293 CELLS

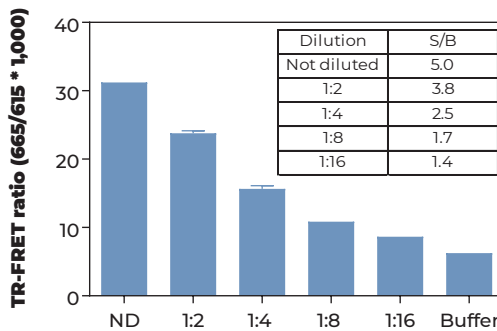


HEK293 cells (50,000 cells/well; in triplicate) were incubated with serial dilutions of the inhibitor BX 795 for 30 min at RT. Cells were then stimulated with 300 nM of Anisomycin for 60 min at RT. Data show that treatment of HEK293 cells with BX 795 inhibits phosphorylation of JNK1/2/3 at T183/Y185 by Anisomycin, but does not affect the levels of total JNK1/2/3.

## HEK293 CONTROL LYSATE TITRATION (QC TEST) PHOSPHO-JNK1/2/3 (T183/Y185)



## HEK293 CONTROL LYSATE TITRATION (QC TEST) TOTAL JNK1/2/3



Quality Control: the Phospho-JNK1/2/3 (T183/Y185) + Total JNK1/2/3 assay kit is routinely tested against Anisomycin-treated HEK293 lysates. HEK293 cells were cultured in a T175 flask to 90% confluence and stimulated with 2  $\mu$ M of Anisomycin 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-JNK1/2/3 (T183/Y185) and total JNK1/2/3. Data show a linear relationship between lysate dilutions and TR-FRET ratio values.



bioauxilium

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