

# Technical Data Sheet

Caution: Research use only.

## Alpha SureFire® Ultra™ Multiplex

### Terbium SureFire Ultra p-JNK 1/2/3 (Thr183/Tyr185) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

**Products: TBSU-PJNK-A500, TBSU-PJNK-A10K, TBSU-PJNK-A50K,  
TBSU-PJNK-A-HV**

#### **Kit-Specificity Information**

This assay kit contains antibodies which recognize the phospho-Thr183/Tyr185 epitope, and a distal epitope, on JNK 1, JNK2 and JNK3. The proteins detected by this kit correspond to GenBank Accessions NP\_620637 (JNK1), NP\_620707 (JNK2) and NP\_620446 (JNK3). JNK1 is also known as SAPK1 and MAPK 8, JNK2 is also known as SAPK, p54aSAPK and MAPK9, and JNK3 is also known as p54bSAPK and MAPK 10. These antibodies recognize JNK 1/2/3 of human, mouse and rat origin. Other species should be tested on a case-by-case basis.

The antibodies in this kit are identical to those in the corresponding ALSU-PJNK-A kits, other than the CaptSure2 tag.

#### **Control Lysate Information**

Positive Control Lysate: HEK293 cells were cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with anisomycin (1µg/mL) for 30min. Cells were then lysed in 10mL of 1X SureFire Ultra Lysis buffer.

#### **Representative data**

Control lysate samples measured in the terbium channel using an Envision plate reader.

