

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ p-JNK 1/2/3 (Thr183/Tyr185) Assay Kit

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

Kit-Specificity Information

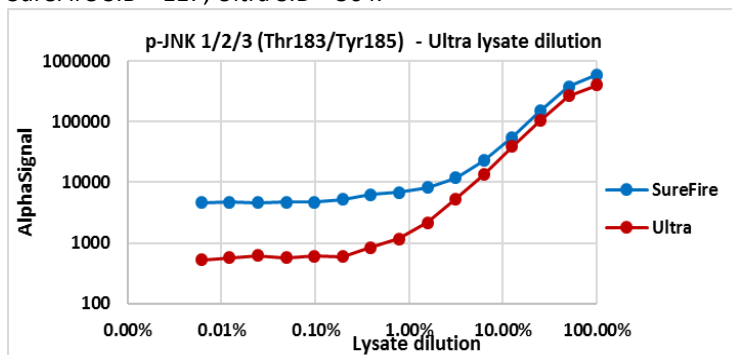
This assay kit contains antibodies which recognize the phospho-Thr183/Tyr185 epitope, and a distal epitope, on JNK 1, JNK2 and JNK3. The proteins detected by this kit correspond to GenBank Accessions NP_620637 (JNK1), NP_620707 (JNK2) and NP_620446 (JNK3). JNK1 is also known as SAPK1 and MAPK 8, JNK2 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.

Control Lysate Information

Positive Control Lysate: Prepared from HEK293 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 1 μ g/mL Anisomycin for 30min and lysed in 10mL of 1X *SureFire Ultra* Lysis buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. Data shows a comparison of AlphaScreen *SureFire* and AlphaLISA *SureFire Ultra* assay performance. Control lysate was diluted in the Lysis buffer specific for the assay type. *SureFire* S:B = 117; *Ultra* S:B = 564.



For more information and detailed protocols, please refer to the generic AlphaLISA *SureFire Ultra* Assay Kits Manual, available from <http://www.perkinelmer.com/category/alpha-surefire-kits>