

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ JNK 1/2 Total Assay Kit

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

Kit-Specificity Information

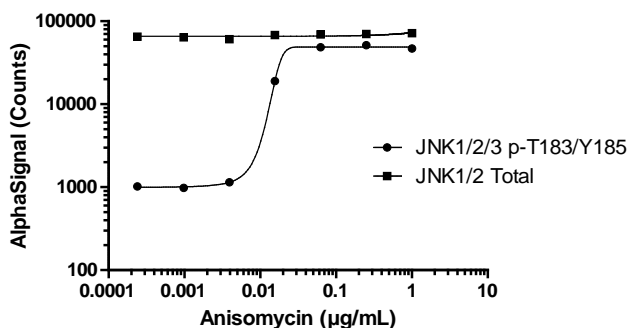
This assay kit contains antibodies which recognize invariant epitopes on JNK 1 and JNK2. The proteins detected by this kit correspond to Entrez Gene ID 5599 and 5601, respectively. JNK1 is also known as PRKM8, SAPK1 and MAPK8 and JNK2 is also known as SAPK, p54aSAPK and MAPK9. These antibodies recognize JNK 1/2 of human 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 µg/mL Anisomycin for 30 min and lysed in 10 mL of 1X *SureFire Ultra* Lysis buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. HEK293 cells were plated overnight at 40K cells/well, starved for 2 hrs, then treated with the indicated concentrations of Anisomycin for 30 mins. Cells were lysed with 100 µL 1X Lysis Buffer and samples assayed separately for total and phospho (T183/Y185) JNK, using respective kits.



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>