

AlphaLISA® SureFire® Ultra™

Human and Mouse p-JNK 1/2/3 (Thr183/Tyr185) Detection Kit

Product number: ALSU-PJNK-A500, ALSU-PJNK-A10K,
 ALSU-PJNK-A50K, ALSU-PJNK-A-HV



Kit specificity:

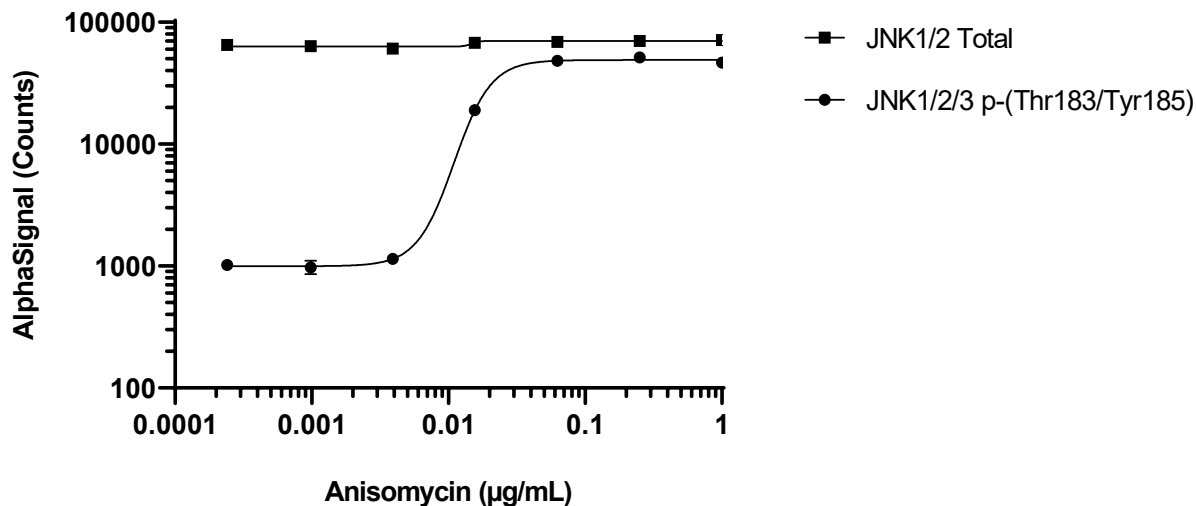
This assay kit contains antibodies which recognize the phospho-Thr183/Tyr185 epitope and a distal epitope, on JNK 1, JNK2 and JNK3. The proteins detected by this kit correspond to UniProt ID P45983 (JNK1), P45984 (JNK2) and P53779 (JNK3). JNK1 is also known as SAPK1c and MAPK8, JNK2 is also known as SAPK1a, and MAPK9, and JNK3 is also known as SAPK1b and MAPK10. These antibodies recognize JNK 1/2/3 of human and mouse origin. Other species should be tested on a case-by-case basis.

Control lysate information:

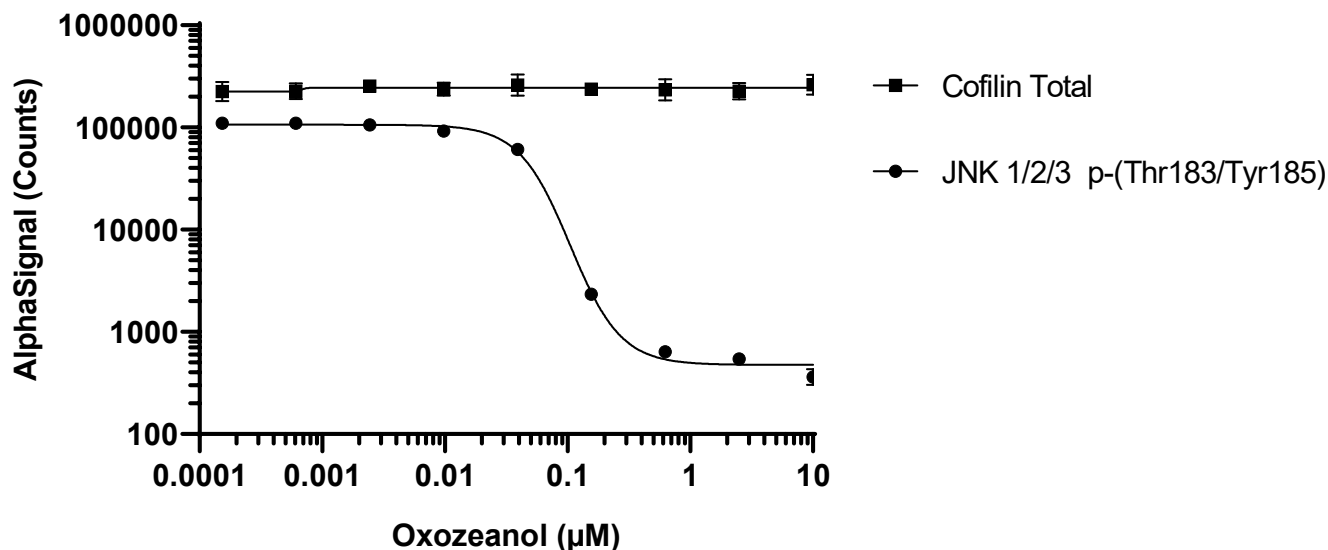
Positive Control Lysate: Prepared from HEK293 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, treated with 1 µg/mL Anisomycin for 30 minutes and lysed with 10 mL of Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. HEK293 cells were plated at 40K cells/well in a 96 well plate and incubated overnight. Cells were serum starved for 2 hours and then treated with Anisomycin at the indicated concentrations for 30 minutes. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Thr183/Tyr185) and Total JNK 1/2, using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint.



Data obtained with a 2-plate, 2-incubation protocol. HEK293 cells were plated at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with Oxozeanol at the indicated concentrations for 2 hours and then 0.75M Sorbitol for 30 minutes. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Thr183/Tyr185) JNK 1/2/3 and Cofilin Total, using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint (JNK p-Thr183/Tyr185) or 400 cells/datapoint (Cofilin Total).



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF"

