

AlphaLISA® SureFire® Ultra™

Human p-MKK4 (Ser257) Detection Kit

Product number: ALSU-PMKK4-A500, ALSU-PMKK4-A10K,
ALSU-PMKK4-A50K, ALSU-PMKK4-A-HV



Kit specificity:

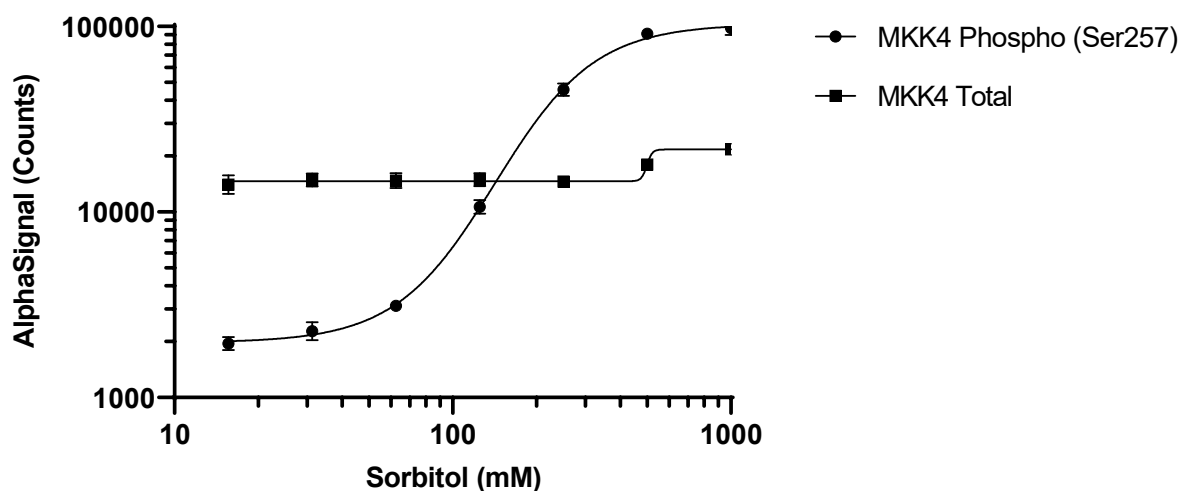
This assay kit contains antibodies which recognize the phospho-Ser257 epitope and a distal epitope, on MKK4. The protein detected by this kit correspond to UniProt ID P45985. MKK4 is also known as MAP2K4, MAPPK 4, MEK4, SEK1 and JNKK1. These antibodies recognize MKK4 of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

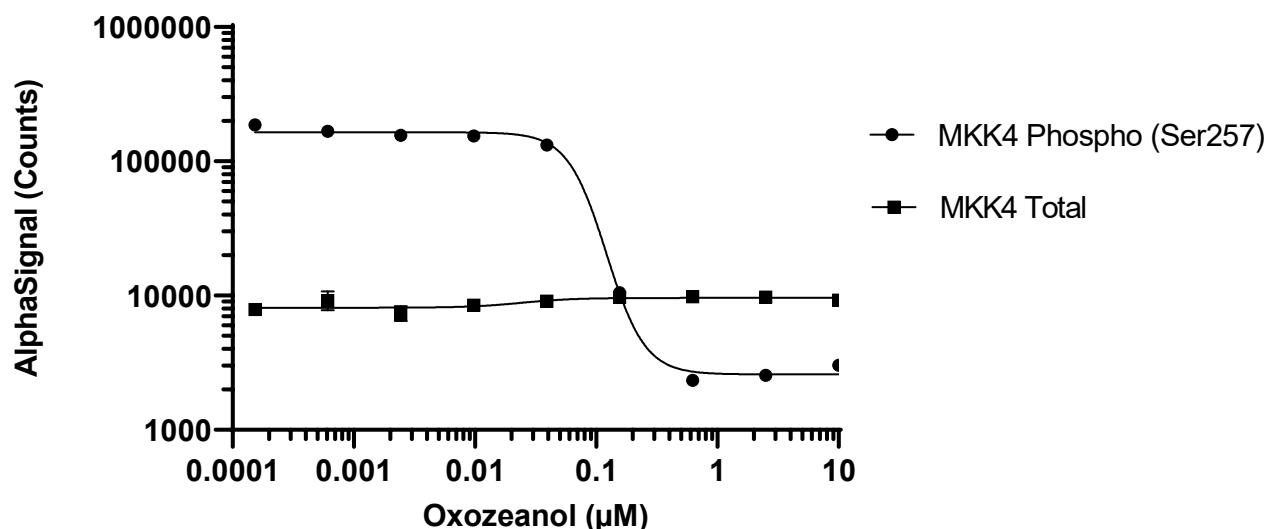
Positive Control Lysate: Prepared from HEK293 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, treated with 1M Sorbitol for 45 minutes and lysed with 4 mL of Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. HEK293 cells were plated at 60K cells/well in a 96 well plate and incubated overnight. Cells were treated with Sorbitol at the indicated concentrations for 45 minutes. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Ser257) and Total MKK4 using respective *SureFire Ultra* kits. Equivalent to approximately 6,000 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 (Ser257) and Total MKK4 using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint.



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"

