

AlphaLISA® SureFire® Ultra™

Human and Mouse MEK2 Total Detection Kit

Product number: ALSU-TMEK2-A500, ALSU-TMEK2-A10K,
 ALSU-TMEK2-A50K, ALSU-TMEK2-A-HV



Kit specificity:

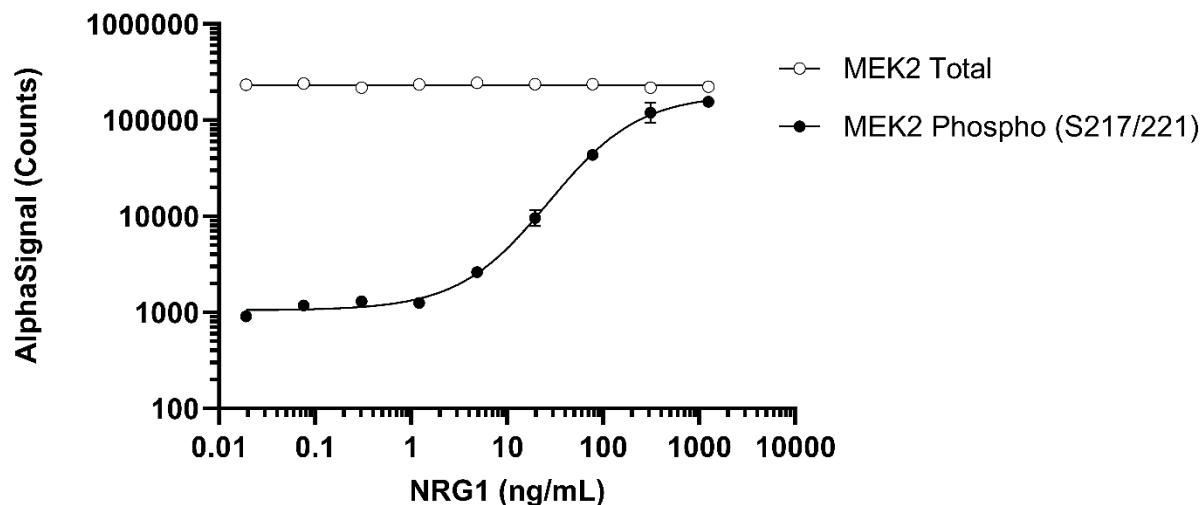
This assay kit contains antibodies which recognize distinct epitopes on MEK2. The protein detected by this kit corresponds to UniProt ID P36507. MEK2 is also known as MAP kinase kinase 2, MAP2K2 and MAPKK2. These antibodies recognize MEK2 of human and mouse origin. Other species should be tested on a case-by-case basis.

Control lysate information:

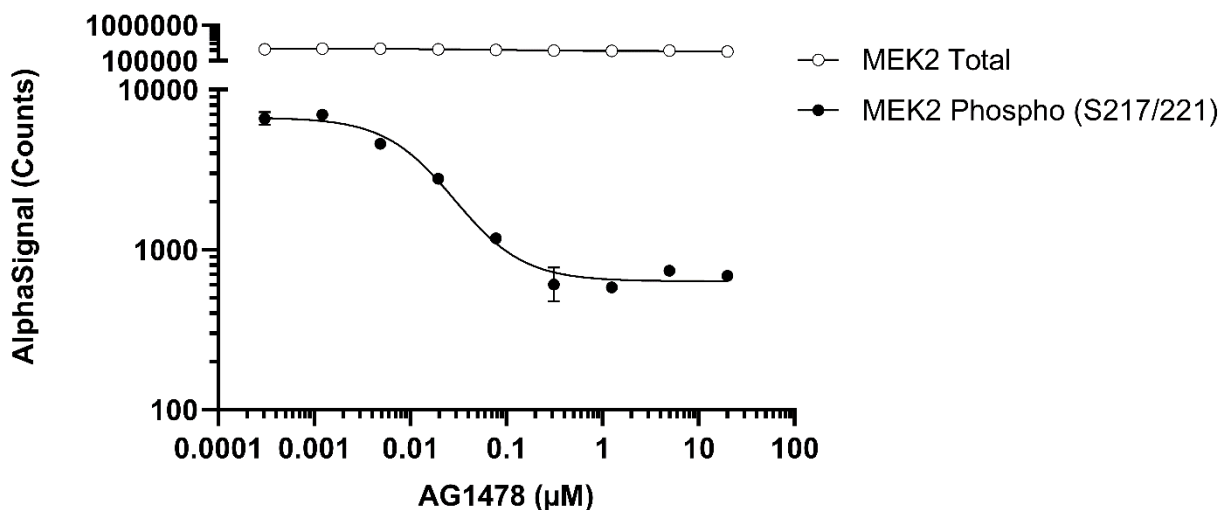
Positive Control Lysate: Prepared from A431 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 200 ng/mL EGF for 10 minutes and lysed with 10 mL of Lysis Buffer.

Representative data

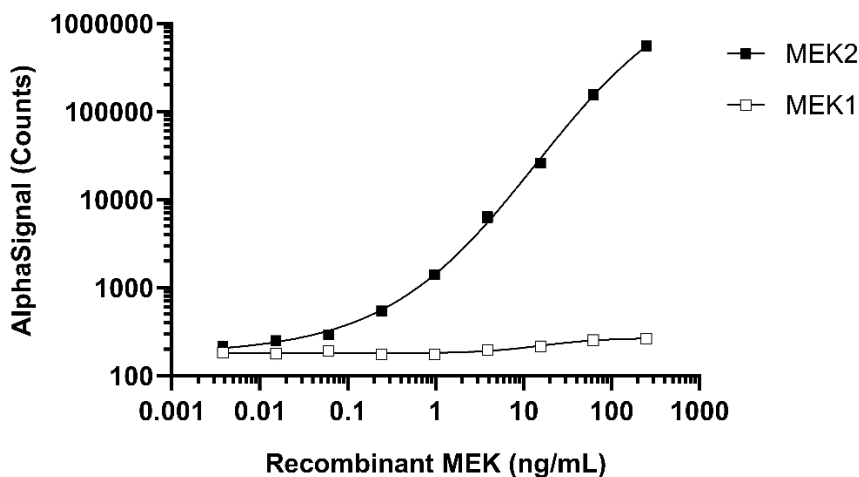
Data obtained with a 2 plate, 2 incubation protocol. MCF7 cells were seeded at 40K cells/well in a 96 well plate and incubated overnight. Cells were starved for 2 hours and then treated with Neuregulin1 (NRG1) at the indicated concentrations for 10 minutes. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Ser117/221) and Total MEK2 using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint.



Data obtained with a 2 plate, 2 incubation protocol. A431 cells were seeded at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with AG1478 at the indicated concentrations for 2 hours and then 1 ng/mL EGF for 30 minutes. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Ser117/221) and Total MEK2 using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint.



MEK1 or MEK2 recombinant proteins were serially diluted in Lysis Buffer and assayed using the Total MEK2 kit. No cross-reactivity against MEK1 recombinant protein was observed.



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"

