

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ p-MEK1 (Ser218/222) Assay Kit

ALSU-PMEK1-A500, ALSU-PMEK1-A10K, ALSU-PMEK1-A50K, ALSU-PMEK1-A-HV

Kit-Specificity

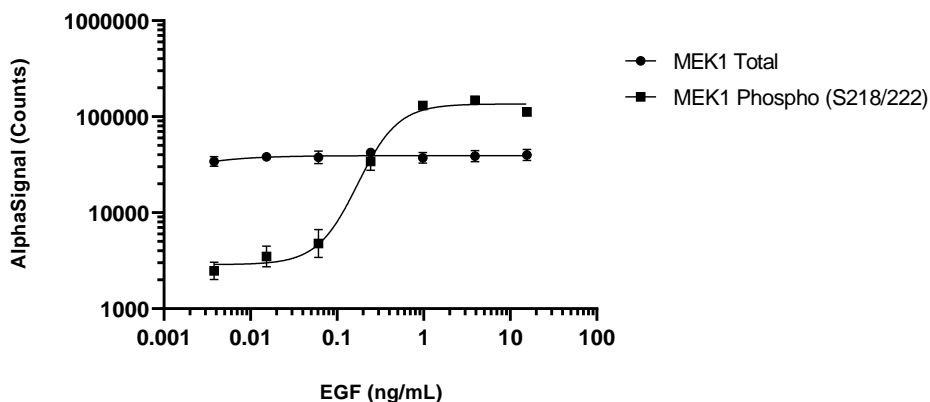
This assay kit contains antibodies which recognizes the phospho-Ser218/222 epitope, and a distal epitope on MEK1. The protein detected by this kit corresponds to UniProt ID Q02750. MEK1 is also known as Dual specificity mitogen-activated protein kinase kinase 1, MKK1 or MAPKK1. These antibodies recognize MEK1 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 in 10 mL of 1X *SureFire Ultra* Lysis Buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. A431 cells were plated at 40K cells per well overnight in a 96 well plate. Cells were then serum starved for 2 hours prior to treatment with EGF at the indicated concentrations for 10 minutes. Cells were lysed with 1x Lysis Buffer and assayed for Total or Phospho (S218/222) MEK1 using respective *SureFire Ultra* kits. Equivalent to approximately 4000 cells/datapoint.



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>