

Alpha SureFire® Ultra™ Multiplex

# Human and Mouse p-MEK1 (Ser218/222) + Total MEK1 Detection Kit

**Product number:** MPSU-PTMEK1-K500, MPSU-PTMEK1-K10K,  
MPSU-PTMEK1-K50K, MPSU-PTMEK1-K-HV



### Kit specificity:

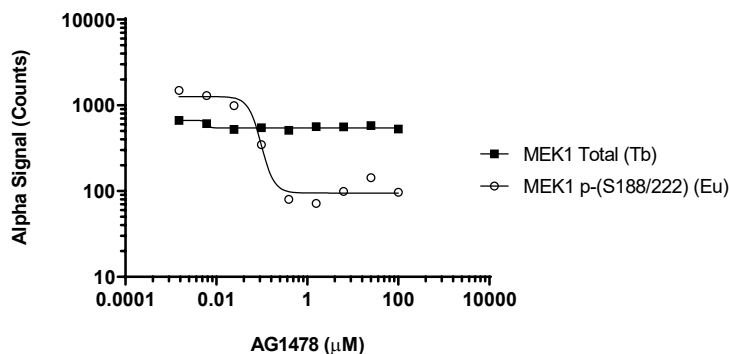
This assay kit contains antibodies which recognize the phospho-Ser218/222 epitope and a distal epitope on MEK1, plus antibodies to detect Total MEK1. The protein detected by this kit corresponds to UniProt ID Q02750. MEK1 is also known as Dual specificity mitogen-activated protein kinase kinase 1, ERK activator kinase 1 and MAP2K1. 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 A-431 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, treated with 200 ng/mL rh-EGF for 10 minutes and lysed with 10 mL of Lysis Buffer.

### Representative data

Data obtained with a 2-plate, 2-incubation protocol. A-431 cells were seeded at 20K cells/well in a 96 well plate and incubated overnight. Cells were treated with AG1478 at the indicated concentrations for 2 hours, followed by 2 ng/mL rh-EGF for 10 minutes. Cells were lysed with Lysis Buffer and assayed for Phospho (Ser218/222) and Total MEK1 in the same wells of an assay plate using the *SureFire Ultra* Multiplex kit. Equivalent to approximately 1,000 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"