

# Technical Data Sheet

Caution: Research use only.

## Alpha SureFire® Ultra™ Multiplex p-MEK1 (Ser218/222) + Total MEK1 Assay Kit

**Products: MPSU-PTMEK1-K500, MPSU-PTMEK1-K10K, MPSU-PTMEK1-K50K,  
MPSU-PTMEK1-K-HV**

### **Kit-Specificity Information**

This assay kit contains antibodies which recognize the phospho-Ser218/222 epitope, and a distal epitope, on MEK1, plus antibodies to measure the total levels of MEK1. The protein detected by this kit corresponds to GenBank Accession NP\_002746. MEK1 is also known as MKK1, MAPKK1, PRKMK1 and MAP2K1. These antibodies recognize MEK1 of human and mouse origin. Other species should be tested on a case-by-case basis.

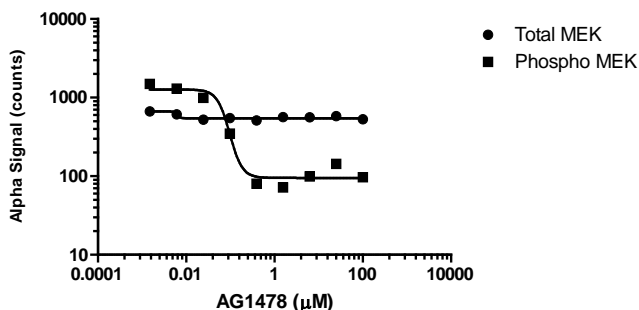
### **Control Lysate Information**

Positive Control Lysate: Prepared from A431 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 200ng/mL EGF for 10min and lysed in 10mL of 1X *SureFire Ultra* Lysis buffer.

### **Representative data**

Data obtained with a 2-plate, 2-incubation protocol. Confluent A431 cells were treated with AG1478 at the indicated concentrations for 2 hours, then stimulated with EGF for 10min. Cells were then lysed with 1X Lysis buffer and assayed for both phospho and total MEK1 as detailed in the MPSU manual.

#### **AG1478 + EGF Dose Response in A431 cells - MEK1 MPSU**



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>