

# Technical Data Sheet

Caution: Research use only.

## AlphaLISA® SureFire® Ultra™ MEK1 Total Assay Kit

**Products: ALSU-TMEK1-A500, ALSU-TMEK1-A10K, ALSU-TMEK1-A50K,  
ALSU-TMEK1-A-HV**

### **Kit-Specificity Information**

This assay kit contains 2 antibodies which recognize invariant epitopes on MEK1. The protein detected by this kit corresponds to GenBank Accession NP\_002746. MEK1 is also known as MKK1, MAPKK1, PRKMK1 and MAP2K1.

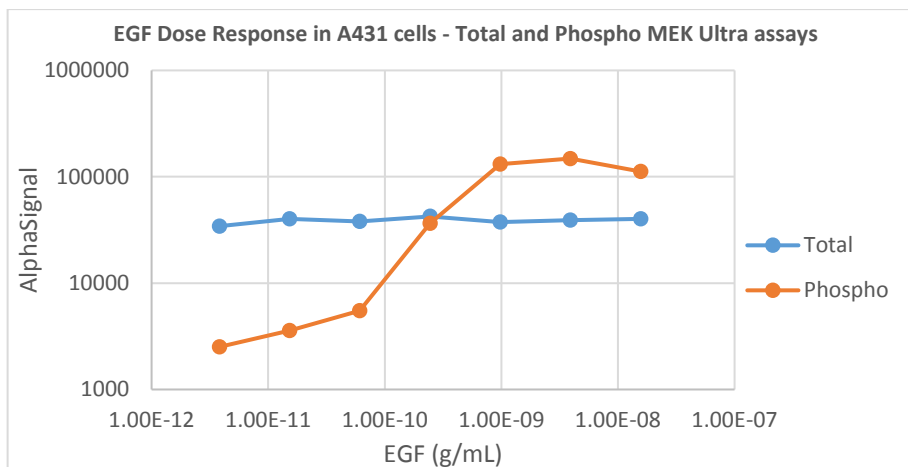
These antibodies recognize MEK1 of human and mouse origin. Other species should be tested on a case-by-case basis.

### **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.



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>