

Technical Data Sheet

Caution: Research use only.

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

**Products: ALSU-PMEK1-A500, ALSU-PMEK1-A10K, ALSU-PMEK1-A50K,
ALSU-PMEK1-A-HV**

Kit-Specificity Information

This assay kit contains 2 antibodies; an antibody which recognizes the phospho-Ser218/222 epitope, and an antibody which recognizes a distal epitope 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.

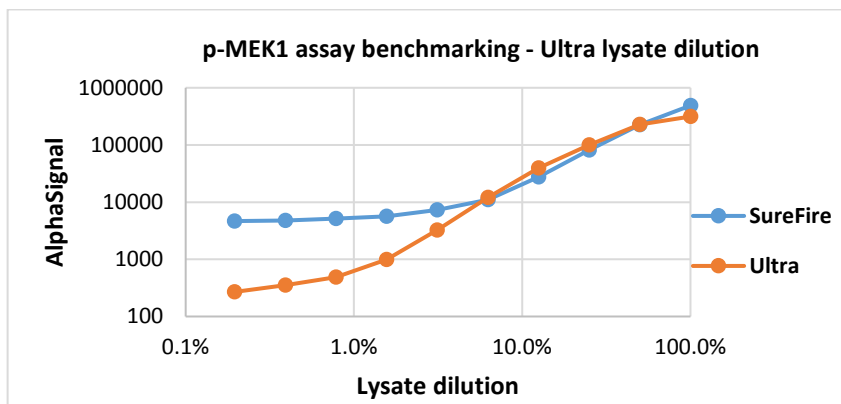
Control Lysate Information

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. Data shows a comparison of AlphaScreen *SureFire* and AlphaLISA *SureFire Ultra* assay performance. Control lysate was diluted in the Lysis buffer specific for the assay type.

S:B *SureFire Ultra* = 1021; S:B *SureFire* = 140



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>

TGRSU176.0