

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex

Terbium SureFire Ultra p-MEK1 (Ser218/222) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

**Products: TBSU-PMEK1-A500, TBSU-PMEK1-A10K, TBSU-PMEK1-A50K,
TBSU-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, PRKM1 and MAP2K1. These antibodies recognize MEK1 of human and mouse origin. Other species should be tested on a case-by-case basis. The antibodies in this kit are identical to those in the corresponding ALSU-PMEK1-A kits, other than the CaptSure2 tag.

Control Lysate Information

Positive Control Lysate: Prepared from A431 cells, cultured to confluence in T175 flasks in 10% FBS containing medium for 3 days, then treated with EGF (200ng/mL) for 10min and lysed in 20mL of 1X SureFire Ultra Lysis buffer.

Representative data

Control lysate samples measured in the terbium channel using an Envision plate reader.

