

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex p-ERK1/2 (Thr202/Tyr204) + Total ERK1/2 Assay Kit

**Products: MPSU-PTERK-M500, MPSU-PTERK-M10K, MPSU-PTERK-M10K,
MPSU-PTERK-M-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Thr202/Tyr204 epitope and a distal epitope, on ERK1/2, plus antibodies to measure the total levels of ERK1/2.

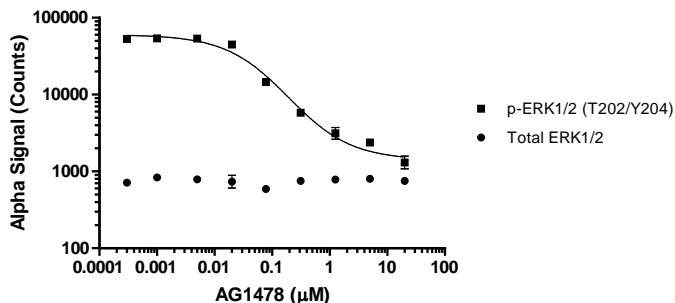
The proteins detected by this kit correspond to Entrez Gene ID 5595,5594. Alternate names include p44 MAPK, MAPK3 (ERK1), p42 MAPK, MAPK1 (ERK2). p44 MAPK and MAPK3 (ERK1), These antibodies recognize ERK1/2 of human, mouse and rat origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

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

Representative data

Data obtained with a 2-plate, 2-incubation protocol. HeLa cells were plated for overnight at 40 K cells/well, then inhibited with the indicated concentrations of AG1478 for 2 hr. Treated with 1 ng/ml EGF for 10 min, then lysed with 100 μ L 1X Lysis Buffer. Samples were assayed for both phospho and total ERK1/2, as detailed in the MPSU manual.



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>