

# Technical Data Sheet

Caution: Research use only.

## AlphaLISA® SureFire® Ultra™ ERK 1/2 Total Assay Kit

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

### **Kit-Specificity Information**

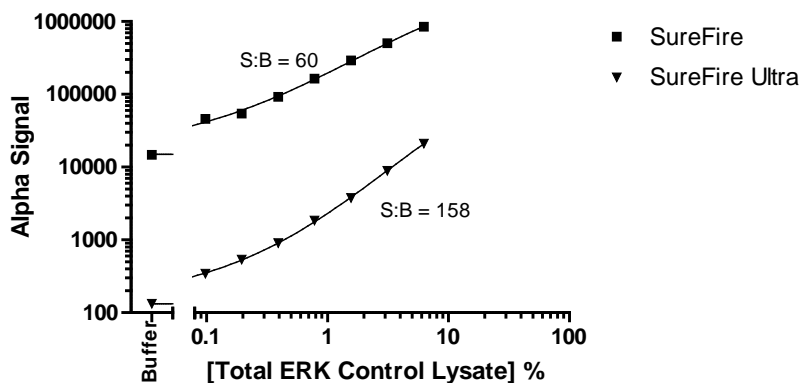
This assay kit contains 2 antibodies which recognize two distinct epitopes on ERK1 and ERK2. The proteins detected by this kit correspond to GenBank Accessions NP 002737 (ERK1) and NP 620407 (ERK2). Alternate Names include p42 MAPK, MAPK3 (ERK1), p42 MAPK, MAPK1 (ERK2). These antibodies recognize ERK 1/2 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 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.



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>