

# Technical Data Sheet

Caution: Research use only.

## AlphaLISA® SureFire® Ultra™ p-AMPK $\alpha$ 1/2 (Thr172) Assay Kit

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

### **Kit-Specificity Information**

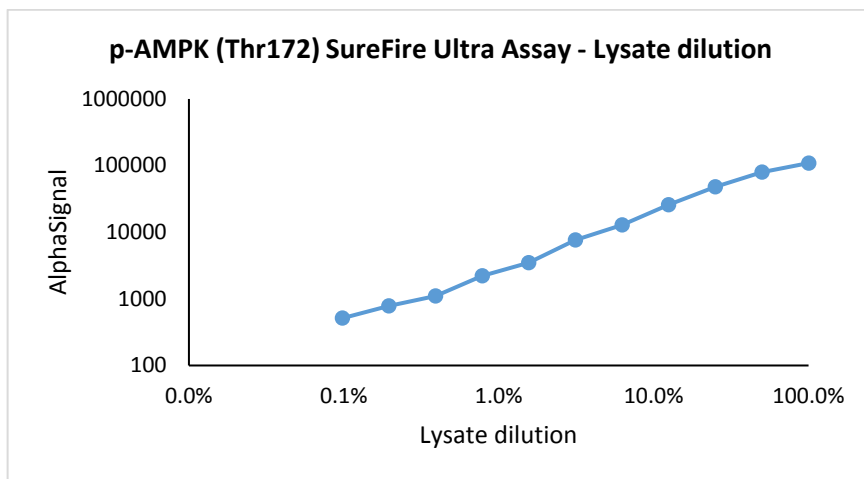
This assay kit contains antibodies which recognize the phospho-Thr172 epitope, and a distal invariant epitope, on p-AMPK $\alpha$ . The assay recognizes both AMPK $\alpha$ 1 and AMPK $\alpha$ 2, corresponding to Entrez-Gene ID's 5562 and 5563, respectively. These antibodies recognize AMPK $\alpha$  of human, mouse and rat origin. Other species should be tested on a case-by-case basis.

### **Control Lysate Information**

Positive Control Lysate: Prepared from HEK293 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 11mM H2O2 for 10min and lysed in 5ml 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>