

# Technical Data Sheet

Caution: Research use only.

## AlphaLISA SureFire® Ultra™

### AMPK $\alpha$ 1/2 Total Assay Kit

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

#### Kit-Specificity Information

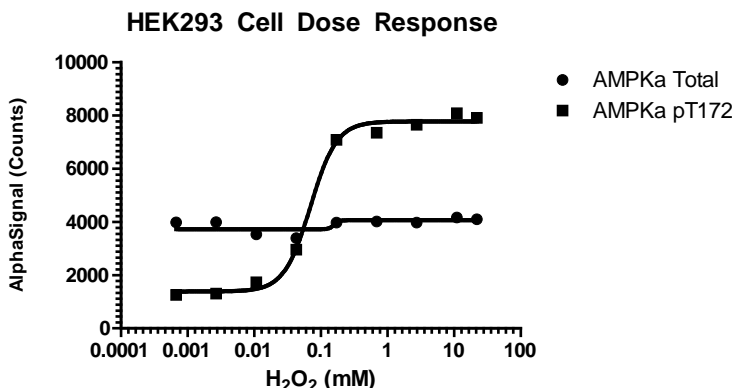
This assay kit contains antibodies which recognize two invariant epitopes on AMPK $\alpha$ . The assay recognizes both AMPK $\alpha$ 1 and AMPK $\alpha$ 2, corresponding to Entrez-Gene ID's 5562 and 5563, respectively. These antibodies recognize AMPK $\alpha$  of human, mouse and rat origin. Other species 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 H<sub>2</sub>O<sub>2</sub> for 10min and lysed in 5ml 1X SureFire Ultra Lysis Buffer.

#### Representative data

Data obtained with a 2-plate, 2-incubation protocol. HEK-293 cells were seeded at 20K cells per well and grown for 2 days in a 96 well plate. Cells were then treated for 10 minutes with peroxide at the indicated concentrations. Cells were lysed with 100 $\mu$ L of 1X Lysis buffer and assayed separately for phospho and total AMPK $\alpha$  with respective SureFire Ultra assay kits.



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>