

# Technical Data Sheet

Caution: Research use only.

## AlphaLISA® SureFire® Ultra™ p-Akt 1/2/3 (Ser473) Assay Kit

**Products: ALSU-PAKT-B500, ALSU-PAKT-B10K, ALSU-PAKT-B50K,  
ALSU-PAKT-B-HV**

### Kit-Specificity Information

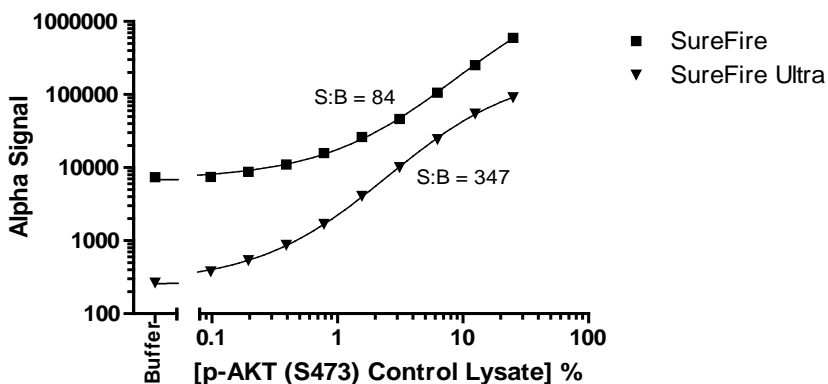
This assay kit contains antibodies which recognize the phospho-Ser473 epitope, and a distal epitope, common to Akt 1, Akt 2 and Akt 3. The proteins detected by this kit correspond to GenBank Accessions NP\_001014431, NP\_001617, and NP\_859029. Akt proteins are also known as PKB or RAC. These antibodies recognize Akt 1/2/3 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 20% FBS for 15min and lysed in 4mL 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>