

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ p-Akt 1 (Ser473) Assay Kit

**Products: ALSU-PAKT1-C500, ALSU-PAKT1-C10K, ALSU-PAKT1-C50K,
ALSU-PAKT1-C-HV**

Kit-Specificity Information

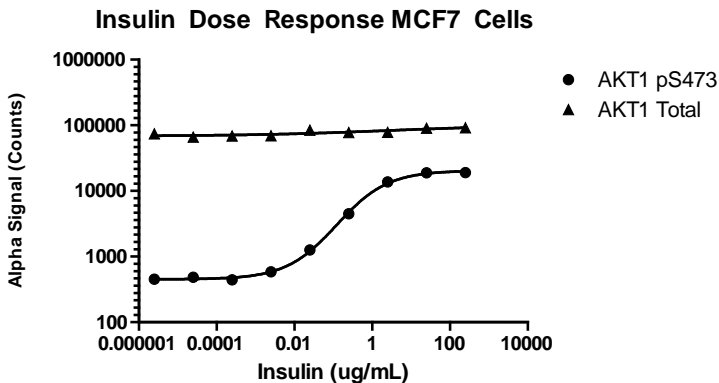
This assay kit contains antibodies which recognize the phospho-Ser473 epitope, and a distal epitope on Akt 1. The kit is not reactive to Akt 2 or Akt 3. The protein detected by this kit correspond to GenBank Accession NP_001014431. Akt proteins are also known as PKB or RAC. These antibodies recognize 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. Confluent MCF-7 cells were serum starved for 4 hours, then stimulated with insulin at the indicated concentrations for 30min. Cells were then lysed with 1X Lysis buffer and samples separately assayed for p-AKT1 (Ser473) and AKT Total, using 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>