

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ Akt1 Total Assay Kit

**Products: ALSU-TAKT1-B500, ALSU-TAKT1-B10K, ALSU-TAKT1-B50K,
ALSU-TAKT1-B-HV**

Kit-Specificity Information

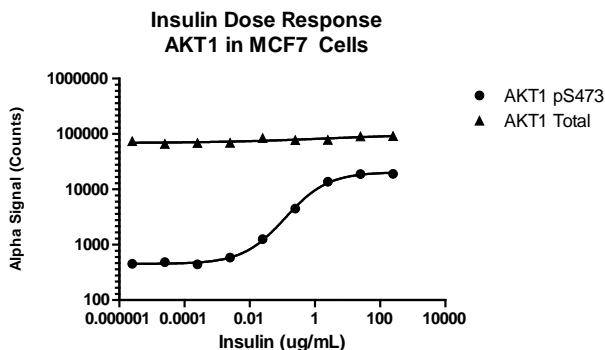
This assay kit contains antibodies which recognize two distinct epitopes on Akt 1. The protein detected by this kit corresponds to GenBank Accession NP_001014431. Also known as PKB, RAC, PRKBA, PKB-ALPHA, RAC-ALPHA. These antibodies recognize Akt 1 of human, mouse and rat origin. Other species should 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 MCF7 cells in a 96 well plate were serum starved for 4 hours, then treated with insulin at the indicated concentrations for 30mins, and cells then lysed with 1X Lysis buffer. Lysates were assayed for total or phospho Akt1 (Ser473) 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>