

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ AKT 1/2/3 Total Assay Kit

Products: ALSU-TAKT-B500, ALSU-TAKT-B10K, ALSU-TAKT-B50K,
ALSU-TAKT-B-HV

Kit-Specificity Information

This assay kit contains antibodies which recognize invariant epitopes on AKT 1/2/3. 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 and mouse origin. Other species should be tested on a case-by-case basis.

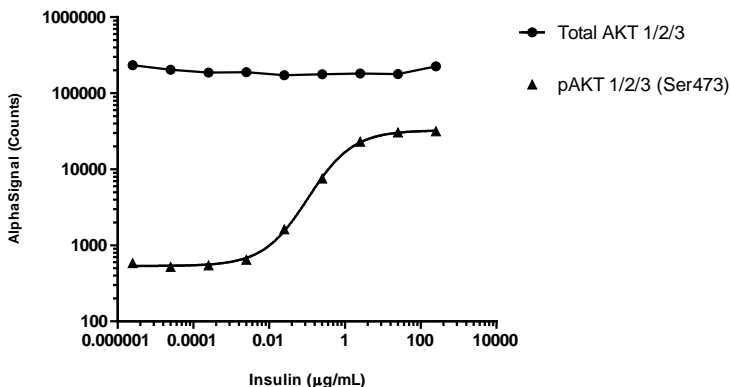
Control Lysate Information

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. MCF-7 cells were plated at 200K cells per mL, incubated overnight, starved for 2 hours, then stimulated with insulin for 30 minutes at the indicated concentrations. Cells were lysed with 100µL of 1X Lysis buffer and assayed.

Insulin Dose Response - MCF7 cells



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>