

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex p-Akt1/2/3 (Ser473) + Total Akt1 Assay Kit

**Products: MPSU-PTAKT-M500, MPSU-PTAKT-M10K, MPSU-PTAKT-M50K,
MPSU-PTAKT-M-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Ser473 epitope, and a distal epitope, on Akt 1/2/3 for phospho Akt analysis, plus antibodies to measure the total levels of Akt 1. The proteins detected by this kit correspond to GenBank Accessions NP_001014431, NP_001617, and NP_859029.

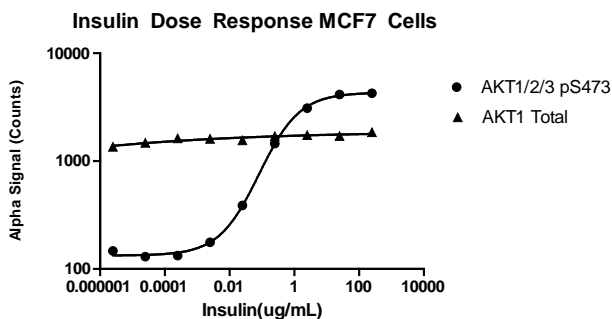
These antibodies recognize Akt of human, mouse and rat origin. Other species should be tested on a case-by-case basis.

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. MCF-7 cells were plated overnight at 20K cells per well in a 96 well plate, serum starved for 2 hours, then treated with insulin for 30 minutes. Cells were lysed with 100µL of 1X Lysis buffer and samples were assayed for both phospho Akt1/2/3 (Ser473) and total Akt1 as detailed in the MPSU manual.



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>