

AlphaLISA® SureFire® Ultra™

Human and Mouse p-AKT 1/2/3 (Thr450) Detection Kit

Product number: ALSU-PAKT-C500, ALSU-PAKT-C10K,
ALSU-PAKT-C50K, ALSU-PAKT-C-HV



Kit specificity:

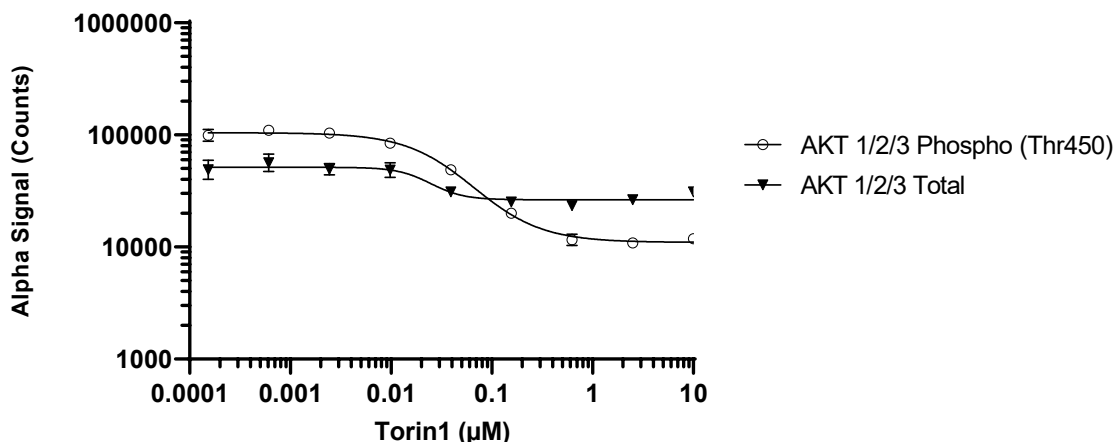
This assay kit contains antibodies which recognize the phospho-Thr450 epitope and a distal epitope on AKT. The protein detected by this kit corresponds to UniProt ID P31749. AKT is also known as PKB or RAC. These antibodies recognize AKT of human and mouse origin. Other species should be tested on a case-by-case basis.

Control lysate information:

Positive Control Lysate: Prepared from MCF7 cells, cultured to confluence in T175 flasks in 10% FBS containing media and lysed with 10 mL of Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. PC3 cells were plated at 20K cells/well in a 96 well plate and incubated over two nights. Cells were treated with Torin 1 at the indicated concentrations for 24 hours. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Thr450) and Total AKT using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint.



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF"