

AlphaLISA™ SureFire® Biotin Free

# Human AKT1/2/3 Total Detection Kit

**Product number:** ASBF-TAKT-A500, ASBF-TAKT-A10K,  
ASBF-TAKT-A50K, ASBF-TAKT-A-HV



**Kit specificity:**

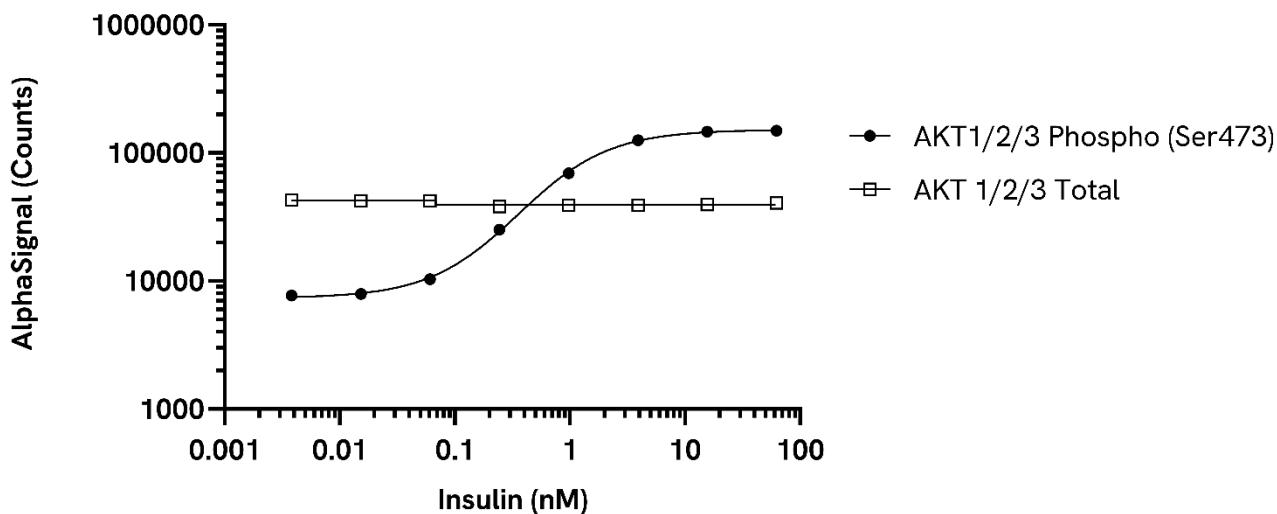
This assay kit contains antibodies which recognize distinct epitopes common to AKT1, AKT2 and AKT3. The proteins detected by this kit corresponds to UniProt ID P31749 (AKT1), P31751 (AKT2) and Q9Y243 (AKT3). AKT proteins are also known as PKB or RAC. These antibodies recognize AKT1/2/3 of human origin. Other species should be tested on a case-by-case basis.

**Control lysate information:**

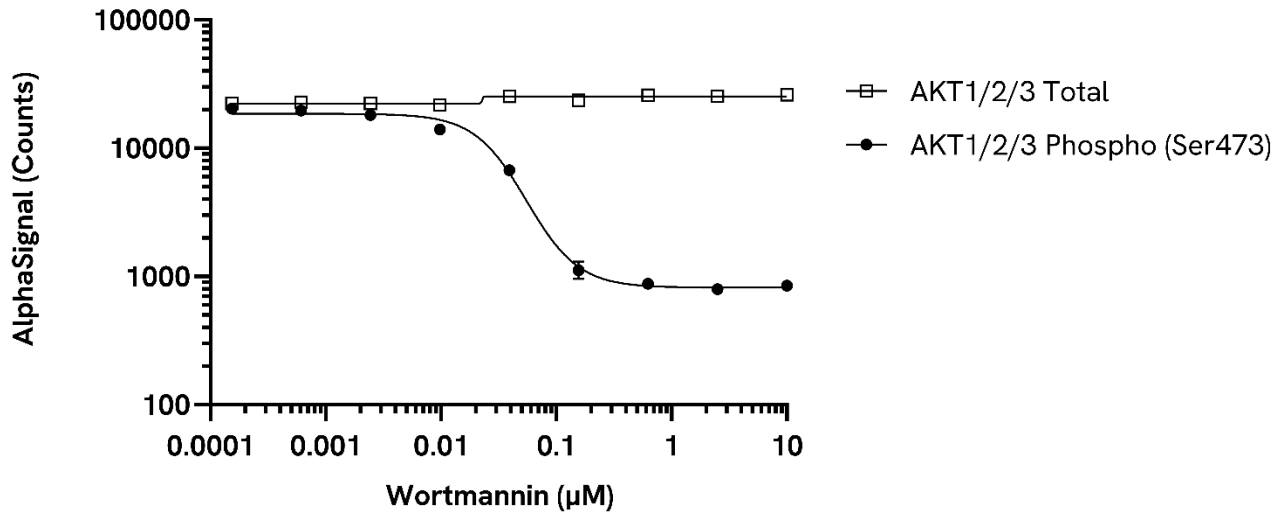
Positive Control Lysate: Prepared from HEK293 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 20% FBS for 15 minutes and lysed with 4 mL of Lysis Buffer.

**Representative data:**

Data obtained with a 2-plate, 2-incubation protocol. THP-1 cells were seeded at 400K cells/well in a 96-well plate and starved in serum free RPMI 1640 medium for 2 hours. Cells were treated with Insulin at the indicated concentrations for 5 minutes. Cells were lysed with the addition of 5X Lysis Buffer and assayed separately for Phospho (Ser473) and Total AKT1/2/3 using respective SureFire Biotin Free kits. Equivalent to approximately 16,000 cells/datapoint.



Data obtained with a 2-plate, 2-incubation protocol. PC3 cells were seeded at 40K cells/well in a 96-well plate in complete Ham's F-12K (Kaighn's) medium and incubated overnight. Cells were treated with Wortmannin at the indicated concentrations in serum free medium for 2 hours. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Ser473) and Total AKT1/2/3 using respective *SureFire* Biotin Free 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"

