

AlphaLISA® SureFire® Ultra™

Human and Mouse RAPTOR Total Detection Kit

Product number: ALSU-TRAPT-A500, ALSU-TRAPT-A10K,
ALSU-TRAPT-A50K, ALSU-TRAPT-A-HV



Kit specificity:

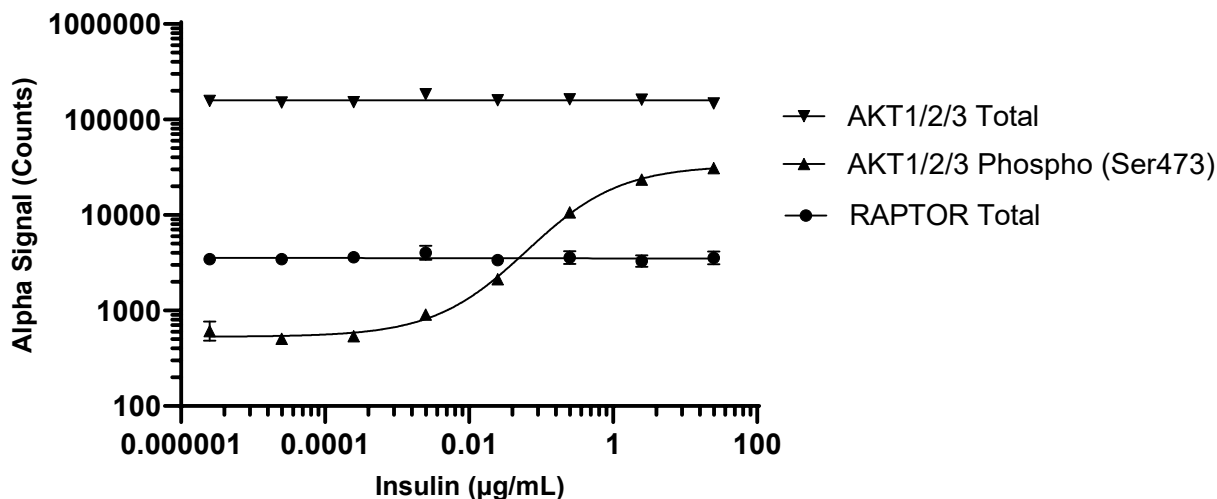
This assay kit contains antibodies which recognize distinct epitopes on RAPTOR. The protein detected by this kit corresponds to UniProt ID Q8N122. RAPTOR is also known as Regulatory-associated protein of mTOR and p150 target of rapamycin (TOR)-scaffold protein. These antibodies recognize RAPTOR of human and mouse origin. Other species should be tested on a case-by-case basis.

Control lysate information:

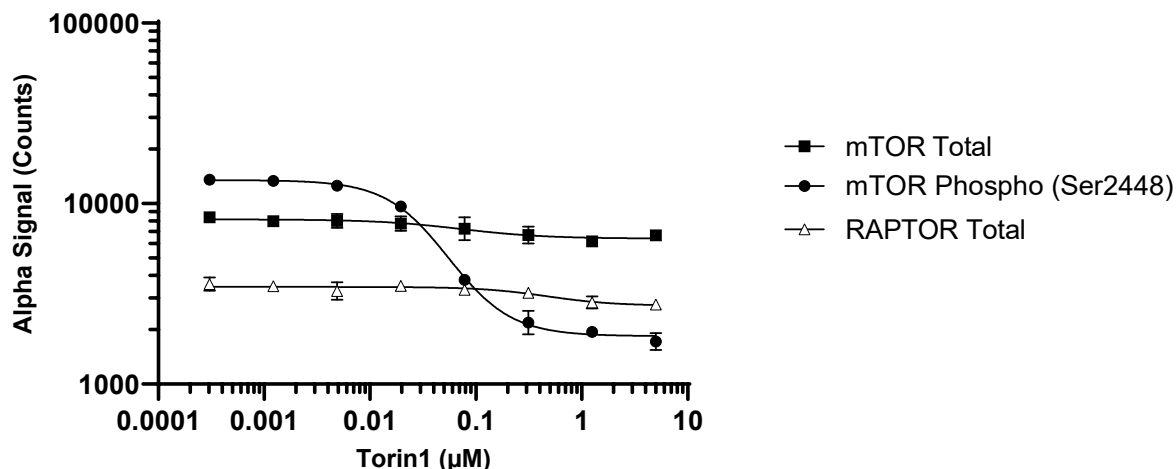
Positive Control Lysate: Prepared from PANC-1 cells, cultured to confluence in T175 flasks in 10% FBS containing media and lysed with 4 mL of Lysis Buffer.

Representative data

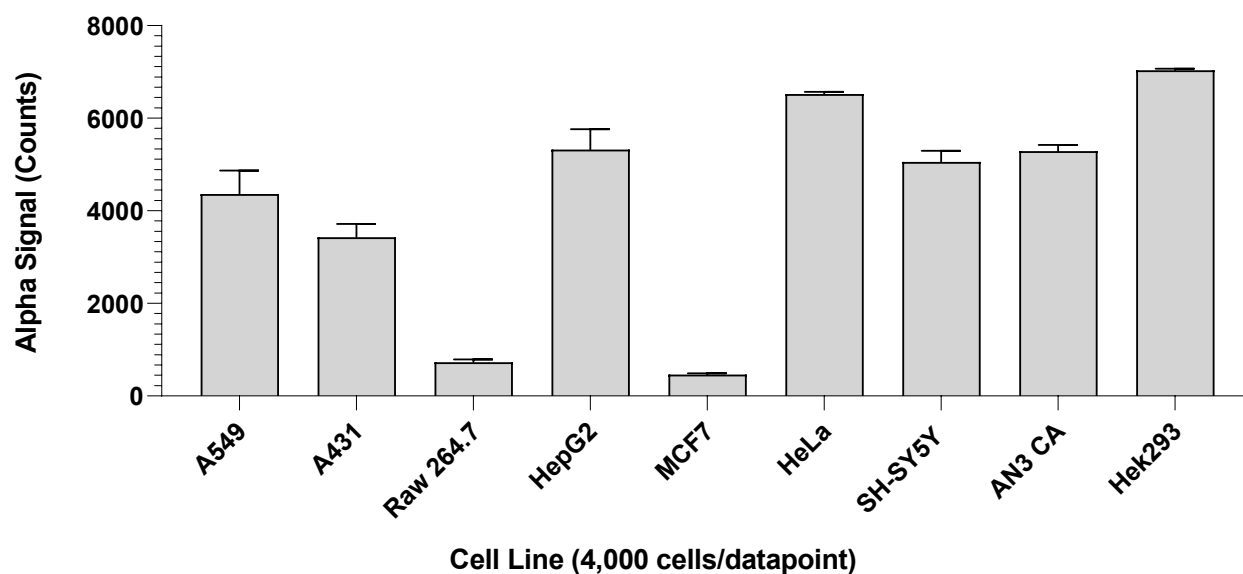
Data obtained with a 2-plate, 2-incubation protocol. MCF7 cells were plated at 40K cells/well in a 96 well plate and incubated overnight. Cells were serum starved for 3 hours prior to treatment with Insulin at the indicated concentrations for 30 minutes. Cells were lysed with Lysis Buffer and assayed separately for various targets using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint.



Data obtained with a 2-plate, 2-incubation protocol. PANC-1 cells were plated at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with a dose range of Torin1 for 18 hours. Cells were lysed with Lysis Buffer and assayed separately for various targets using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint.



Data obtained for Raptor Total from screening various lysates prepared from cells lysed in Lysis Buffer. Approximate number of cells/datapoint is indicated for the various cell lines.



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"

