

AlphaLISA® SureFire® Ultra™

Human and Mouse p-p70S6K (Thr389) Detection Kit

Product number: ALSU-PP70-A500, ALSU-PP70-A10K,
 ALSU-PP70-A50K, ALSU-PP70-A-HV



Kit specificity:

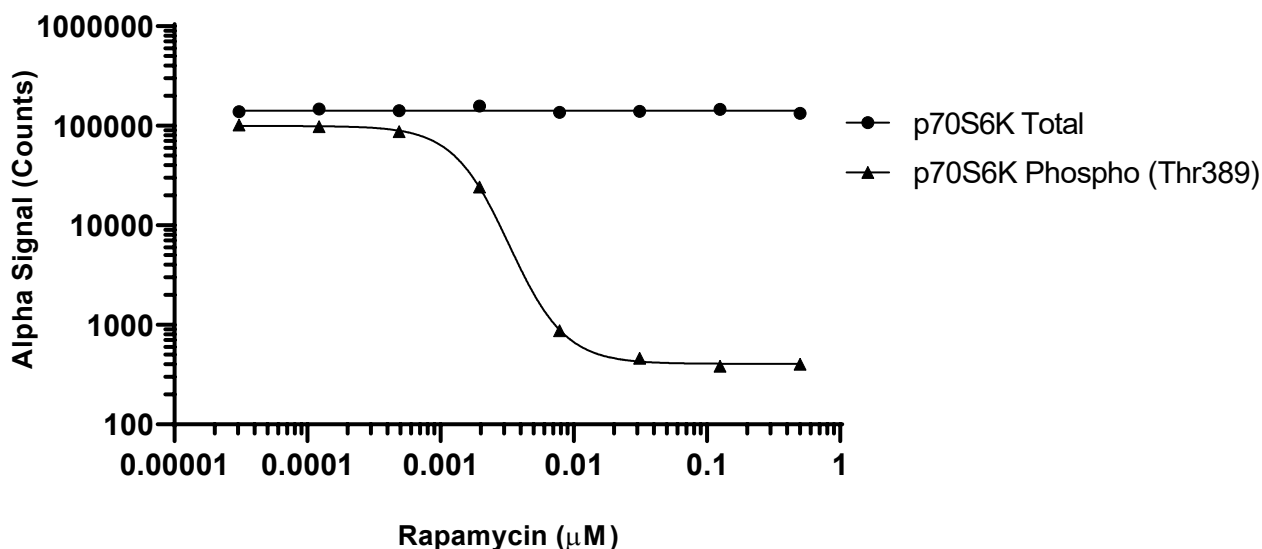
This assay kit contains antibodies which recognize the phospho-Thr389 epitope and a distal epitope on p70S6K. The protein detected by this kit corresponds to UniProt ID P23443. Also known as Ribosomal protein S6 Kinase beta-1 and S6K1. These antibodies recognize p70S6K of human and mouse origin. Other species should be tested on a case-by-case basis.

Control lysate information:

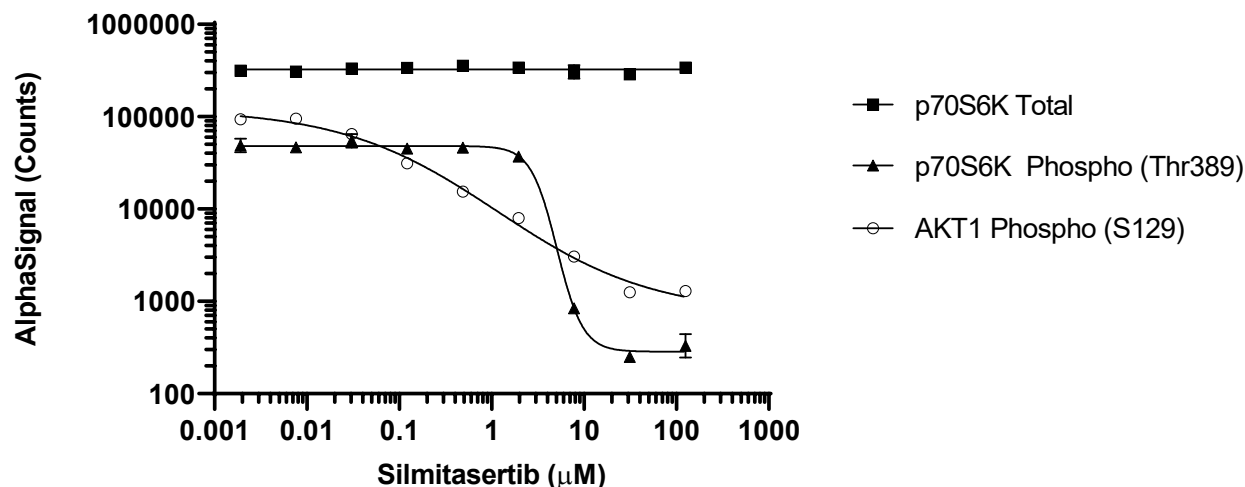
Positive Control Lysate: Prepared from MCF7 cells, cultured to confluence in T175 flasks in 10% FBS containing media, then treated with 500 µg/mL Insulin for 30 minutes and lysed with 20 mL of Lysis Buffer.

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 treated with Rapamycin at the indicated concentrations for 2 hours before the addition of 10 µg/mL Insulin for 30 minutes. Cells were lysed with Lysis Buffer and assayed for Phospho (Thr389) and Total p70S6K using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint.



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 treated with Silmitasertib (CK2 inhibitor) in serum free medium at the indicated concentrations for 24 hours. Cells were lysed with Lysis Buffer and assayed for Phospho (Thr389), Total p70S6K and Phospho (Ser129) AKT1 using respective *SureFire Ultra* kits. Equivalent to approximately 2,000 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"

