

Alpha SureFire® Ultra™ Multiplex

Human and Mouse p-Ribosomal Protein S6 (Ser240/244) + Total Ribosomal Protein S6 Detection Kit

Product number: MPSU-PTRPS6-A500, MPSU-PTRPS6-A10K,
MPSU-PTRPS6-A50K, MPSU-PTRPS6-A-HV



Kit specificity:

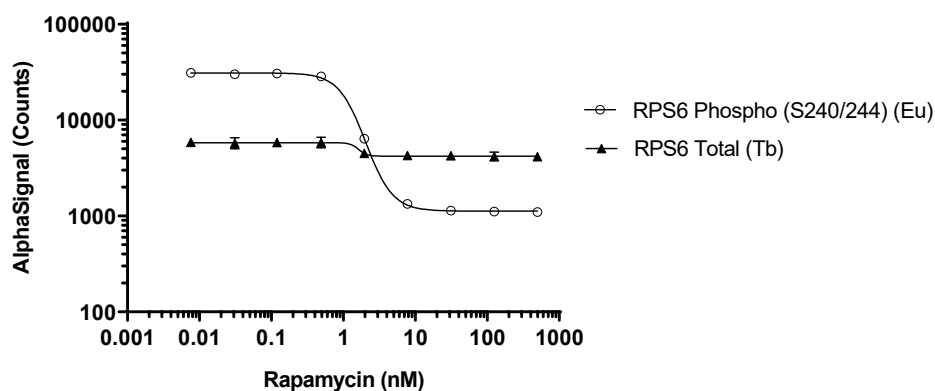
This assay kit contains antibodies which recognize the phospho-RPS6 (Ser240/244) epitope and a distal epitope on RPS6, plus antibodies to detect Total RPS6. The protein detected by this kit corresponds to UniProt ID P62753. These antibodies recognize RPS6 of human and mouse origin. Other species should 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 medium, starved for 2 hours, then treated with 500 µg/mL Insulin for 60 minutes and lysed in 50 mL of Lysis buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. MCF7 cells were seeded at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with Rapamycin at the indicated concentrations for 2 hours and then 2.5 µg/mL Insulin for a further 30 minutes. Cells were lysed with Lysis buffer and assayed for Phospho (Ser240/244) and Total RPS6 in the same wells of an assay plate using the *SureFire Ultra* Multiplex kit. 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"