

AlphaLISA® SureFire® Ultra™ Terbium

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

Product number: TBSU-PS6R-A500, TBSU-PS6R-A10K
 TBSU-PS6R-A50K, TBSU-PS6R-A-HV



Kit Specificity

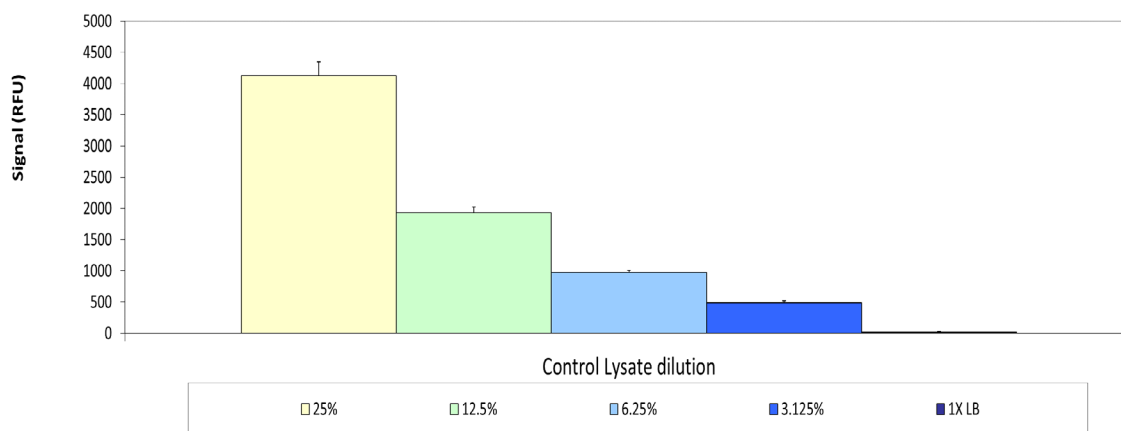
This assay kit contains antibodies which recognize the phospho-Ser240/244 epitope and a distal epitope on Ribosomal Protein S6. The protein detected in this kit corresponds to UniProt ID P62753. RPS6 is also known as Small ribosomal subunit protein eS6 and 40S ribosomal protein S6. These antibodies recognize 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, serum starved for 2 hours then treated with 500 µg/mL Insulin for 1 hour and lysed with 50mL of Lysis Buffer.

Representative data

Data obtained from the positive control lysate diluted in Lysis Buffer. Samples were assayed for Phospho Ribosomal Protein S6 (Ser240/244) using the Terbium SureFire Ultra kit.



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"