

# Technical Data Sheet

## AlphaLISA® SureFire® Ultra™ p70S6K Total Assay Kit

**ALSU-TP70-A500, ALSU-TP70-A10K, ALSU-TP70-A50K, ALSU-TP70-A-HV**

### **Kit-Specificity**

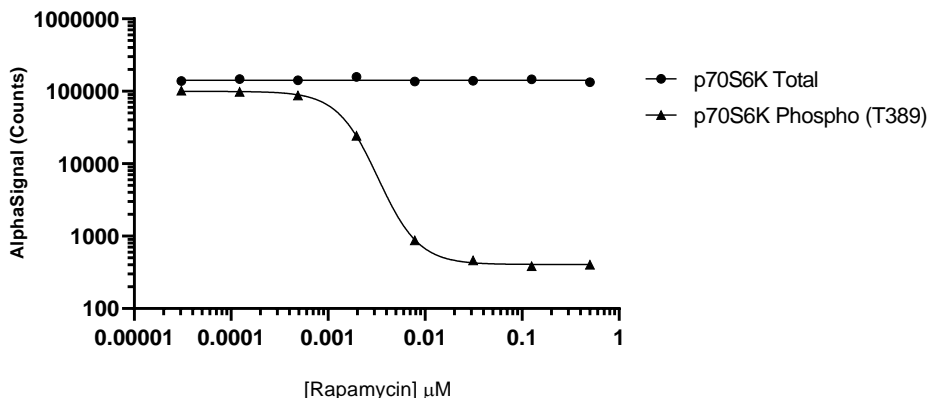
This assay kit contains antibodies which recognize two invariant epitopes on p70S6K. The protein detected by this kit corresponds to UniProt ID P23443. Also known as Ribosomal protein S6 Kinase beta-1, S6K1, p70(S6K)-alpha, and RPS6KB1. These antibodies recognize p70S6K of human, mouse and rat origin. 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, then treated with 500 µg/mL insulin for 30 min and lysed in 20 mL of 1X *SureFire Ultra* Lysis Buffer.

### **Representative Data**

Data obtained with a 2-plate, 2-incubation protocol. MCF7 cells were plated @ 40K cells per well in a 96 well plate and incubated overnight. Cells were then treated with Rapamycin at the indicated concentrations for 2 hours before the addition of 10 µg/mL of Insulin for 30 minutes. Cells were lysed with 1x Lysis Buffer and assayed for Total or Phospho (T389) p70S6K with respective *SureFire Ultra* kits. Equivalent to approximately 4000 cells/datapoint.



For more information and detailed protocols, please refer to the generic AlphaLISA *SureFire Ultra* Assay Kits Manual, available from <http://www.perkinelmer.com/category/alpha-surefire-kits>