

# Technical Data Sheet

## AlphaLISA® SureFire® Ultra™ p-p70S6K (Thr389) Assay Kit

ALSU-PP70-A500, ALSU-PP70-A10K, ALSU-PP70-A50K, ALSU-PP70-A-HV

### Kit-Specificity

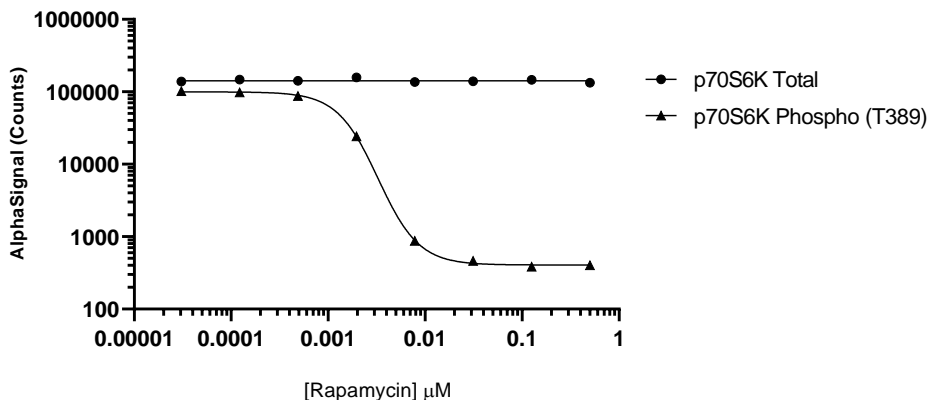
This assay kit contains antibodies which recognize the phospho-Thr389 epitope, and a distal epitope, on p70S6K. The protein detected by this kit corresponds to UniProt ID P23443. Also known as Ribosomal protein S6 Kinase beta-1, S6K1, p70(S6K)-alpha, and RPS6KB1. These antibodies recognize p70S6K of 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 @ at 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>