

Technical Data Sheet

Caution: Research use only.

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

**Products: ALSU-PP70-A500, ALSU-PP70-A10K, ALSU-PP70-A50K,
ALSU-PP70-A-HV**

Kit-Specificity Information

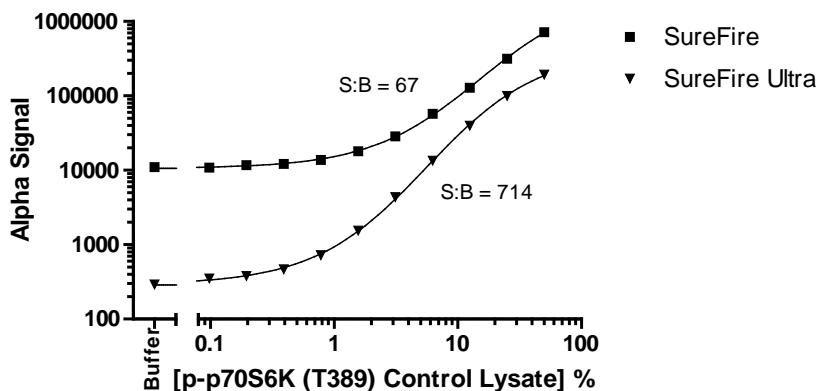
This assay kit contains antibodies which recognize the phospho-Thr389 epitope, and a distal epitope, on p70S6K. The protein detected by this kit corresponds to GenBank Accession NP_003152. Also known as S6K, S6K1, p70-alpha, p70(S6K)-alpha, and RPS6KB1. These antibodies recognize p70S6K of human, mouse 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 500µg/mL Insulin for 30min and lysed in 20mL of 1X *SureFire Ultra* Lysis buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. Data shows a comparison of AlphaScreen *SureFire* and AlphaLISA *SureFire Ultra* assay performance. Control lysate was diluted in the Lysis buffer specific for the assay type.



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>