

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex

Terbium SureFire Ultra p-p70 S6K (Thr389) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

Products: TBSU-PP70-A500, TBSU-PP70-A10K, TBSU-PP70-A50K, TBSU-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.

The antibodies in this kit are identical to those in the corresponding ALSU-PP70-A kits, other than the CaptSure2 tag.

Control Lysate Information

Positive Control Lysate: Prepared from MCF7 cells, cultured to confluence in T175 flasks in 10% FBS containing medium for 3 days, then treated with insulin (500µg/mL) for 30min and lysed in 30mL of 1X SureFire Ultra Lysis buffer.

Representative data

Control lysate samples measured in the terbium channel using an Envision plate reader.

