

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex

Terbium SureFire Ultra p-Ribosomal Protein S6 (Ser240/244) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

**Products: TBSU-PS6R-A500, TBSU-PS6R-A10K, TBSU-PS6R-A50K,
TBSU-PS6R-A-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-RPS6 (Ser240/244) epitope, and a distal epitope, on RPS6. The protein detected by this kit corresponds to GenBank Accession NP_001001. These antibodies recognize RPS6 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-PS6R-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 60min and lysed in 25mL of 1X SureFire Ultra Lysis buffer.

Representative data

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

