

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex

Terbium SureFire Ultra p-mTOR (Ser2448) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

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

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Ser2448 epitope, and a distal epitope, on mammalian target of rapamycin (mTOR). The protein detected by this kit corresponds to GenBank Accession NP_004949. mTOR is also known as FRAP, FRAP1, FRAP2, RAFT1, RAPT1, FLJ44809 and MTOR. These antibodies recognize mTOR 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-PMTOR-A kits, other than the CaptSure2 tag.

Control Lysate Information

Positive Control Lysate: Prepared from Jurkat cells, cultured in a T175 flask to approximately 106 cells/mL in 10% FBS containing medium, then treated with EGF (2µg/mL) for 30min and lysed in 4mL of 1X SureFire Ultra Lysis buffer.

Representative data

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

