

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ p-mTOR (Ser2448) Assay Kit

**Products: ALSU-PMTOR-A500, ALSU-PMTOR-A10K, ALSU-PMTOR-A50K,
ALSU-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.

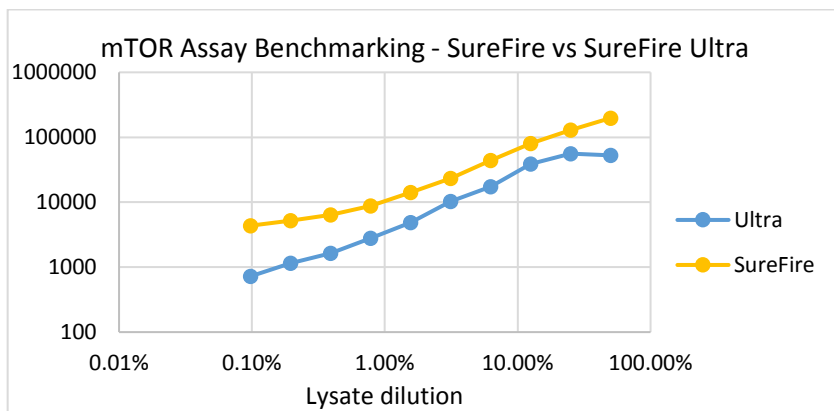
Control Lysate Information

Positive Control Lysate: Prepared from Jurkat cells, cultured in a T175 flask to approximately 1×10^6 cells/mL in 10% FBS containing medium, then treated with $2 \mu\text{g/mL}$ EGF for 30min and lysed at 1×10^7 cells/mL with 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.

S:B *SureFire* = 28; S:B *SureFire Ultra* = 110.



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>