

Technical Data Sheet

Caution: Research use only.

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

**Products: ALSU-PMTOR-C500, ALSU-PMTOR-C10K, ALSU-PMTOR-C50K,
ALSU-PMTOR-C-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 and mouse origin. Other species should be tested on a case-by-case basis.

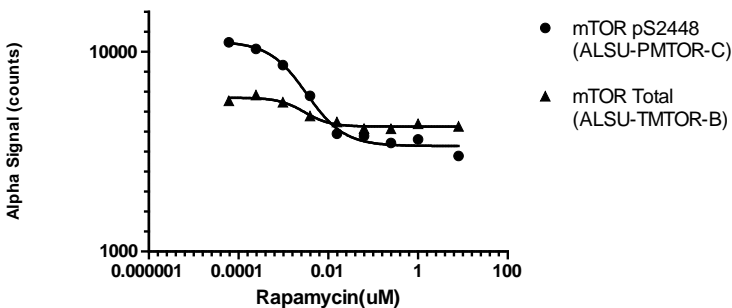
Control Lysate Information

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. HEK cells were plated at 60K cells per well, and treated with Rapamycin for 2 hours at the indicated concentrations, then lysed with 1X Lysis buffer. Levels of mTOR (Ser2448) phosphorylation, and total mTOR were assayed with respective assay kits.

HEK293 Cell Inhibition Dose Response



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>