

# Technical Data Sheet

Caution: Research use only.

## AlphaLISA® SureFire® Ultra™ mTOR Total Assay Kit

**Products: ALSU-TMTOR-A500, ALSU-TMTOR-A10K, ALSU-TMTOR-A50K,  
ALSU-TMTOR-A-HV**

### **Kit-Specificity Information**

This assay kit contains antibodies which recognize two invariant epitopes 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.

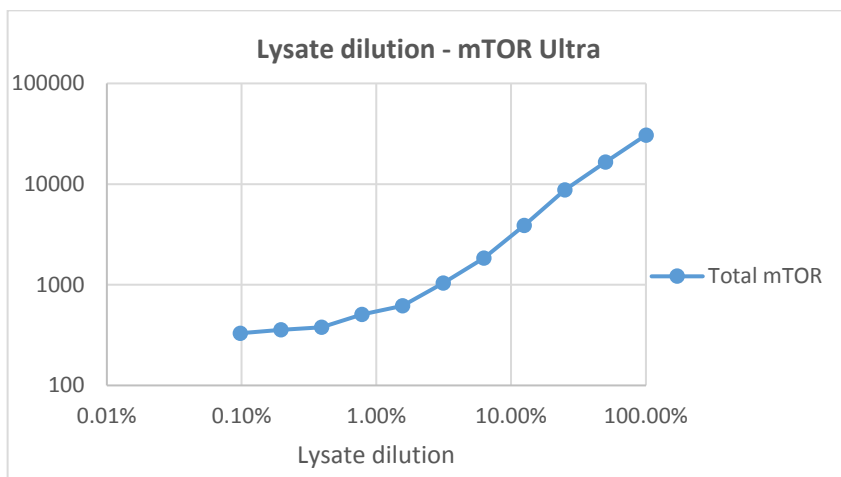
These antibodies recognize mTOR of human, mouse and rat origin. Other species should be tested on a case-by-case basis.

### **Control Lysate Information**

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

### **Representative data**

Data obtained with a 2-plate, 2-incubation protocol.



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-kit>