

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ MLKL Total (Mouse) Assay Kit

ALSU-TMLKL-B500, ALSU TMLKL-B10K, ALSU-TMLKL-B50K, ALSU-TMLKL-B-HV

Kit Specificity

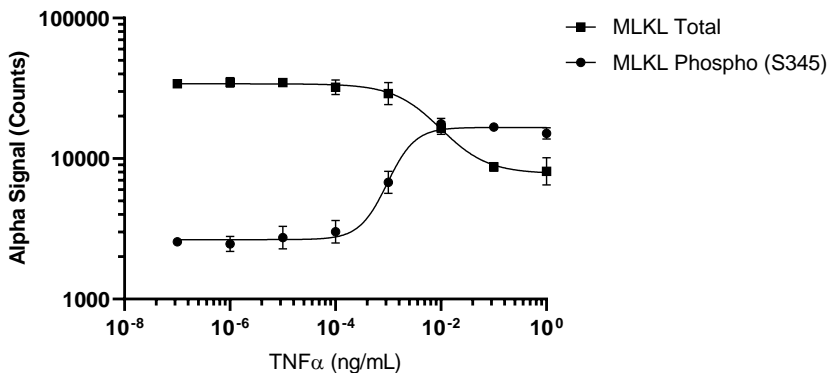
This assay kit contains antibodies which recognize invariant epitopes on MLKL. The protein detected by this kit corresponds to UniProt ID Q8NB16. MLKL is also known as Mixed lineage kinase domain-like protein. These antibodies recognize MLKL of human and mouse origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive control lysate: Prepared from L929 cells, cultured to confluence in T175 flasks in 10% FBS containing medium. Cells were treated with 20 μ M Z-VAD, 100 nM SM-164 and 50 ng/mL TNF α for 4 hours at 37°C. Cells were lysed in 5 mL 1X *SureFire Ultra* Lysis buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. L929 cells were seeded at 4×10^4 cells per well in a 96 well plate and incubated overnight. Cells were treated with 20 μ M Z-VAD, 100 nM SM-164 and TNF α at the indicated concentrations for 4 hours at 37°C. Cells were lysed with 1X *SureFire Ultra* Lysis buffer, and assayed for Phospho (Ser345) or Total MLKL in separate wells of an assay plate with respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint.



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>