

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ p-MLKL (Ser345) (Mouse) Assay Kit

ALSU-PMLKL-A500, ALSU-PMLKL-A10K, ALSU-PMLKL-A50K, ALSU-PMLKL-A-HV

Kit Specificity

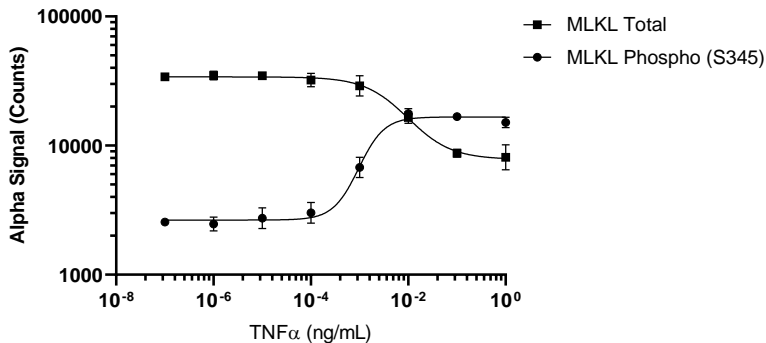
This assay kit contains antibodies which recognize the phospho-Ser345 epitope, and a distal epitope, on MLKL. The protein detected by this kit corresponds to UniProt ID Q8NB16. MLKL is also known as Mixed lineage kinase domain-like protein. These antibodies recognize MLKL of mouse origin. Other species 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>