

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ p38 MAPK Total Assay Kit

ALSU-TP38-B500, ALSU-TP38-B10K, ALSU-TP38-B50K, ALSU-TP38-B-HV

Kit-Specificity

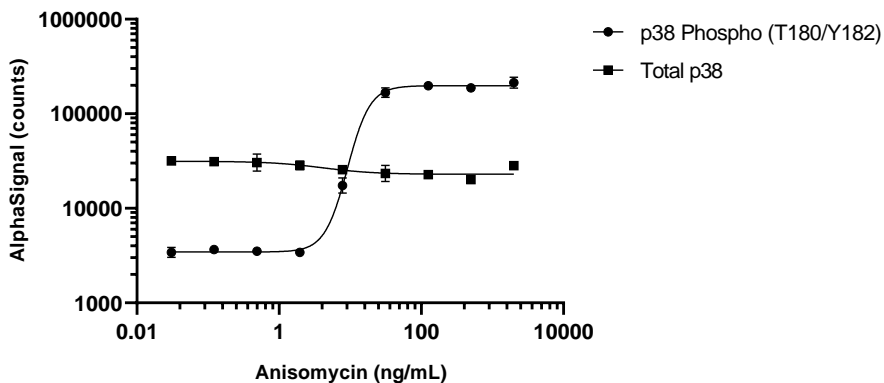
This assay kit contains antibodies which recognize invariant epitopes on p38 MAPK. The protein detected by this kit corresponds to UniProt ID Q16539. P38 MAPK is also known as MAPK14, SAPK2a and MAX-interacting protein 2. These antibodies recognize p38 MAPK of human, mouse, rat origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from HeLa cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 2 µg/mL Anisomycin for 15 min and lysed in 5 mL of 1X *SureFire Ultra* Lysis Buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. HeLa cells were plate at 40K cell per well in a 96 well plate and incubated overnight. Cells were serum starved for 2 hours prior to treatment with Anisomycin at the indicated concentrations for 15 minutes. Cells were lysed with 1x Lysis Buffer and assayed separately for Total or Phospho (T180/Y182) p38 using respective *SureFire Ultra* kits. Equivalent to approximately 4000 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>