

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ p-p38 MAPK (Thr180/Tyr182) Assay Kit

**Products: ALSU-PP38-B500, ALSU-PP38-B10K, ALSU-PP38-B50K,
ALSU-PP38-B-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Thr180/Tyr182 epitope, and a distal epitope, on p38 MAPK. The protein detected by this kit corresponds to GenBank Accession NP_001306. Also known as p38; SAPK2A; p38alpha; MAPK14. These antibodies recognize p38 MAPK of human, mouse, rat origin. Other species should be tested on a case-by-case basis.

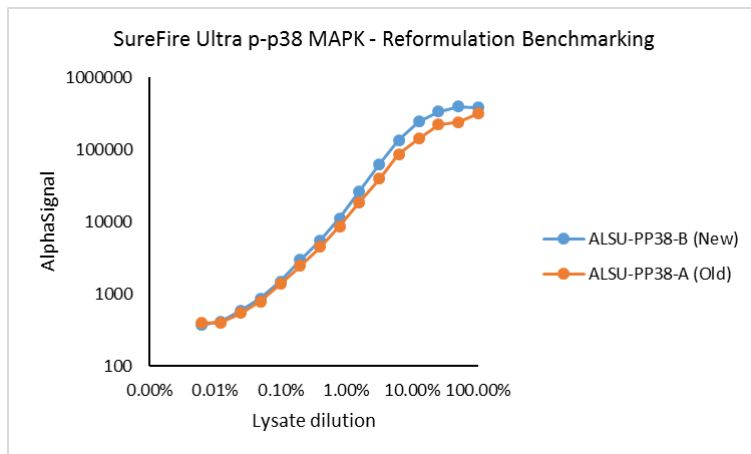
Control Lysate Information

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.

Data is shown for this new formulation kit and the previous old formulation kit.



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>
TGRSU181.0