

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ Human p-HSP27 (Ser15) Detection Kit

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

Kit Specificity

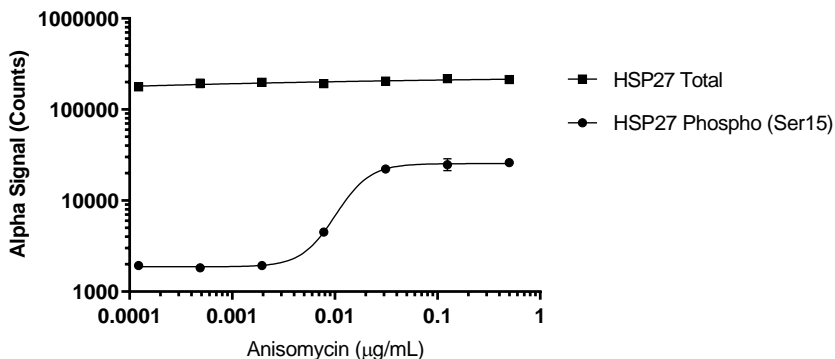
This assay kit contains antibodies which recognize the phospho-Ser15 epitope and a distal epitope on HSP27. The protein detected by this kit corresponds to UniProt ID P04792. HSP27 is also known as heat shock protein family B (small) member 1, HSPB1 and HSP28. These antibodies recognize HSP27 of human origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from A549 cells grown to confluence in T175 flasks in 10% FBS containing medium, then treated with 100 ng/mL Anisomycin for 30 minutes at 37°C. Cells were lysed in 8 mL of *SureFire Ultra* Lysis Buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. A549 cells were plated overnight at 10K cells/well in a 96 well plate, then treated with Anisomycin at the indicated concentrations for 30 minutes. Cells were lysed with *SureFire Ultra* Lysis Buffer and assayed separately for Phospho (Ser15) and Total HSP27, using respective *SureFire Ultra* kits. Equivalent to approximately 1,000 cells/datapoint.



For more information and detailed protocols, please refer to the generic AlphaLISA *SureFire Ultra* Detection Kits Manual, available from <http://www.perkinelmer.com/category/alpha-surefire-kits>