

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ p-STAT1 (Ser727) Assay Kit

ALSU-PST1-B500, ALSU-PST1-B10K, ALSU-PST1-B50K, ALSU-PST1-B-HV

Kit-Specificity

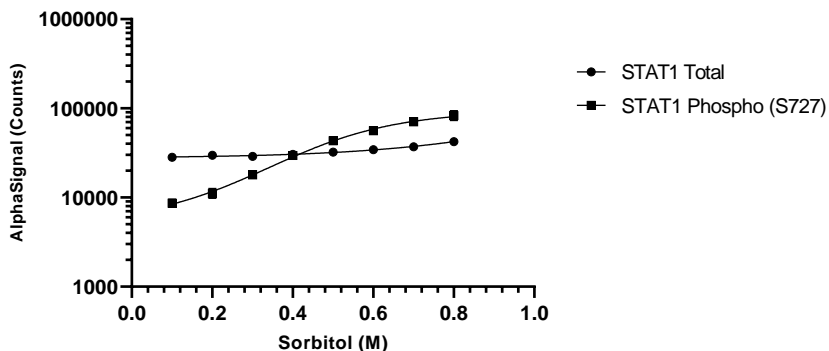
This assay kit contains antibodies which recognize the phospho-Ser727 epitope, and a distal epitope on STAT1. The protein detected by this kit corresponds to UniProt ID P42224. STAT1 is also known as signal transducer and activator of transcription-1. These antibodies recognize STAT1 of human, mouse and 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 0.6 M Sorbitol and 200 ng/mL IFN- γ for 20 minutes each and lysed in 6 mL of 1X *SureFire Ultra* Lysis Buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. HeLa cells were plated at 40K cells per well in a 96 well plate and incubated overnight. Cells were treated with D-Sorbitol at the indicated concentrations for 20 minutes prior to a separate treatment with 200 ng/mL IFN- γ for 20 minutes. Cells were lysed with 1x Lysis Buffer and assayed separately for Total or Phospho (Ser727) Stat1 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>