

Technical Data Sheet

Alpha SureFire® Ultra™ Multiplex p-STAT1 (Ser727) + Total STAT1 Assay Kit

MPSU-PTST1-M500, MPSU-PTST1-M10K, MPSU-PTST1-M50K, MPSU-PTST1-M-HV

Kit-Specificity Information

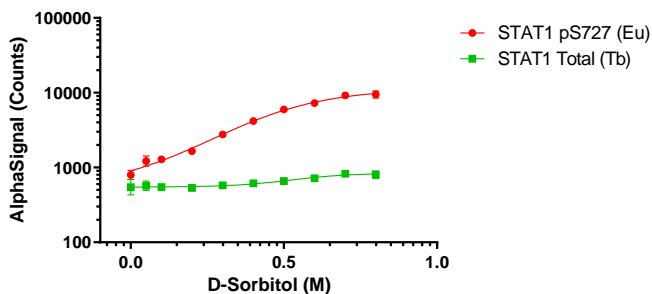
This assay kit contains antibodies which recognize the phospho STAT1 Ser-727 epitope and invariant epitopes on STAT1. The protein detected by this kit corresponds to UniProt ID P42224. STAT1 is also known as signal transducer and activator of transcription-1. These antibodies recognize STAT1 of human 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 consecutively with 0.6 M Sorbitol and 200 ng/mL Interferon-gamma for 20 minutes (independent treatments) 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 seeded overnight at 40K cells/well in a 96 well plate in MEM + 10% FBS. Cells were pre-treated with a dose range of sorbitol for 20 mins prior to a separate 20 min treatment with 200 ng/mL IFN γ . Cells were lysed with 1X *SureFire Ultra* Lysis buffer and assayed for Phospho (Tyr727) and Total STAT1 in the same wells of an assay plate by Alpha *SureFire Ultra* Multiplex. 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>