

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ p-STAT1 (Tyr701) Assay Kit

**Products: ALSU-PST1-A500, ALSU-PST1-A10K, ALSU PST1-A50K,
ALSU-PST1-A-HV**

Kit-Specificity Information

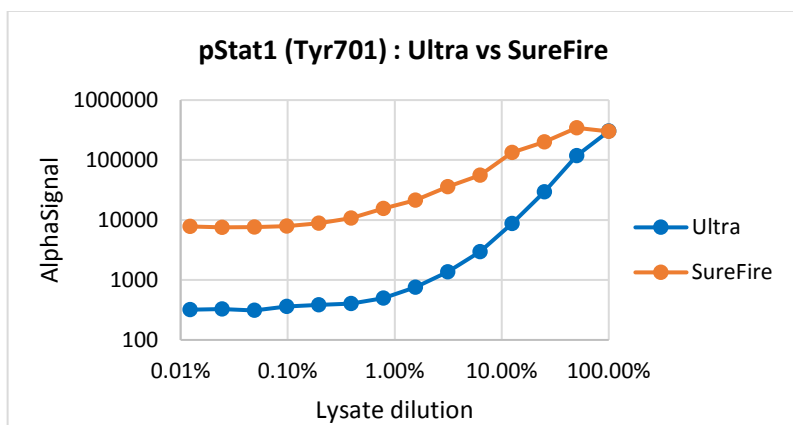
This assay kit contains antibodies which recognize the phospho-Tyr701 epitope, and a distal epitope, on signal transducer and activator of transcription 1 (STAT1). The protein detected by this kit corresponds to GenBank Accession NP_009330. STAT1 is also known as ISGF-3, STAT1 and DKFZp686B04100. These antibodies recognize STAT1 of human and mouse 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 0.6M Sorbitol and 200ng/mL Interferon-gamma for 20min each and lysed in 6mL of 1X *SureFire Ultra* Lysis buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. Data shows a comparison of AlphaScreen *SureFire* and AlphaLISA *SureFire Ultra* assay performance. Control lysate was diluted in the Lysis buffer specific for the assay type.



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>