

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ p-STAT3 (Tyr705) Assay Kit

**Products: ALSU-PST3-A500, ALSU-PST3-A10K, ALSU-PST3-A50K,
ALSU-PST3-A-HV**

Kit-Specificity Information

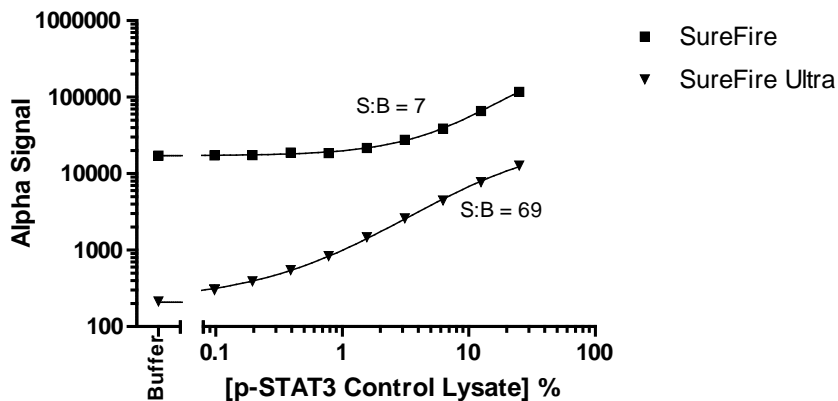
This assay kit contains antibodies which recognize the phospho-Tyr705-epitope, and a distal epitope on STAT3 α and STAT3 β . The proteins detected by this kit correspond to GenBank Accessions NP_644805 and NP_998827. These antibodies recognize STAT3 of human, mouse and rat origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from A431 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 2 μ g/mL EGF for 10min and lysed in 5mL 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>