

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ p-STAT6 (Tyr641) Assay Kit

**Products: ALSU-PST6-A500, ALSU-PST6-A10K, ALSU-PST6-A50K,
ALSU-PST6-A-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Tyr641 epitope, and a distal epitope, on STAT6.

These antibodies recognize STAT6 of human origin. It is predicted to react also with mouse. Other species should be tested on a case-by-case basis

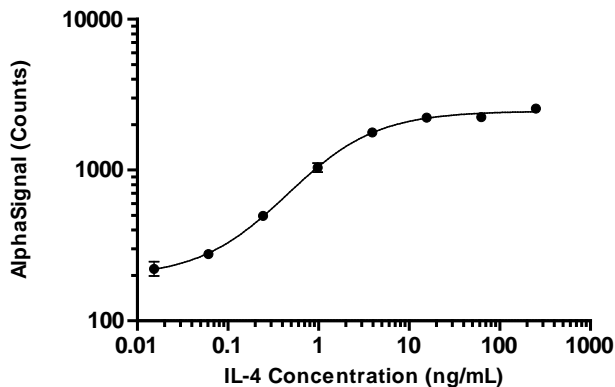
Control Lysate Information

Positive Control Lysate: Prepared from HeLa cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 20 ng/mL rhIL-4 for 20 mins and lysed in 4mL of 1X *SureFire Ultra* Lysis buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. Graph below shows data from HeLa cells plated overnight at 200K cells/mL, treated with IL-4 at concentrations shown for 20 min and lysed in 1X *SureFire Ultra* Lysis buffer.

pSTAT6 (Tyr 641) SF Ultra - IL-4 Dose Response



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>