

# Technical Data Sheet

Caution: Research use only.

## AlphaLISA® SureFire® Ultra™ p-STAT4 (Tyr693) Assay Kit

**Products: ALSU-PST4-A500, ALSU-PST4-A10K, ALSU-PST4-A50K,  
ALSU-PST4-A-HV**

### **Kit-Specificity Information**

This assay kit contains antibodies which recognize the phospho-Tyr693 epitope, and a distal epitope, on STAT4. The protein detected by this kit corresponds to Entrez-Gene Id [6775](#).

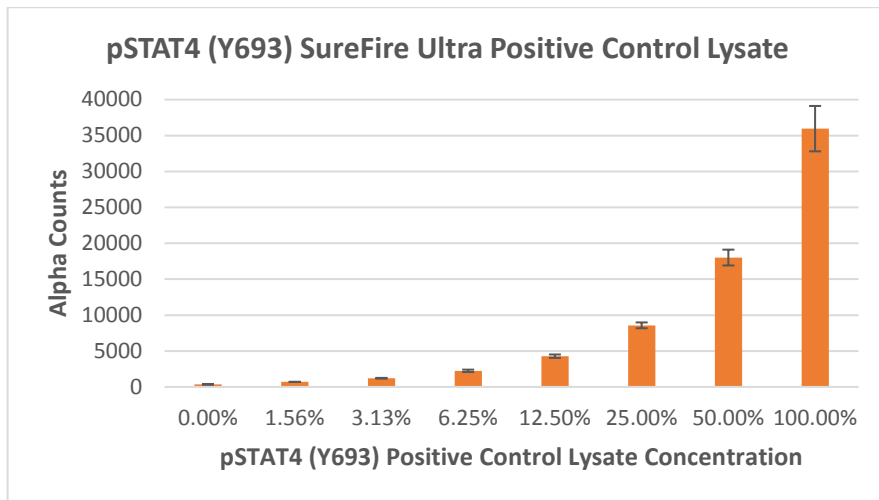
These antibodies recognize STAT4 of human origin. It is predicted to react also with mouse. Other species should be tested on a case-by-case basis.

### **Control Lysate Information**

Positive Control Lysate: Prepared from NK92 cells, cultured to  $0.2 \times 10^6$  cells/mL in T175 flasks in 12.5% FBS containing medium, IL-2 starved, then treated with 200ng/mL IL-12 for 30min and lysed at  $3.2 \times 10^6$  cells/mL with 1X *SureFire Ultra* Lysis buffer.

### **Representative data**

Data obtained with a 2-plate, 2-incubation protocol.



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>