

# Technical Data Sheet

Caution: Research use only.

## Alpha SureFire® Ultra™ Multiplex p-STAT3 (Tyr705) + Total STAT3 Assay Kit

**Products: MPSU-PTST3-K500, MPSU-PTST3-K10K, MPSU-PTST3-K50K,  
MPSU-PTST3-K-HV**

### **Kit-Specificity Information**

This assay kit contains antibodies which recognize the phospho-Tyr705 epitope, and a distal epitope, on STAT3 for phospho STAT3 analysis, plus antibodies to measure the total levels of STAT3.

The proteins detected by this kit correspond to Entre Gene ID 6774. These antibodies recognize STAT3 of human, mouse and rat origin. Other species should be tested on a case-by-case basis.

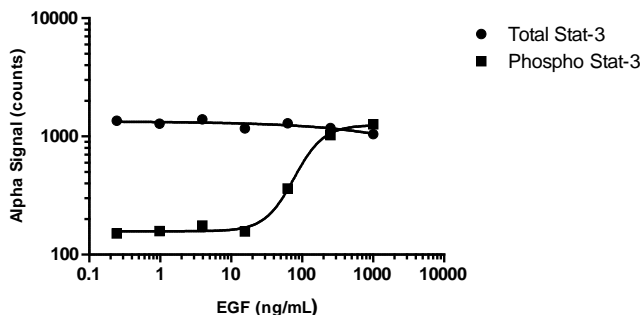
### **Control Lysate Information**

Positive Control Lysate: Prepared from A431 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 2µg/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. Confluent A431 cells were stimulated with EGF at the indicated concentrations for 10min, then lysed with 1X Lysis buffer and assayed for both phospho and total STAT3 as detailed in the MPSU manual.

#### **EGF Dose Response in A431 cells - Stat3 MPSU**



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>