

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ p-STAT5 (Tyr694/699) Assay Kit

ALSU-PST5-B500, ALSU-PST5-B10K, ALSU-PST5-B50K, ALSU-PST5-B-HV

Kit-Specificity

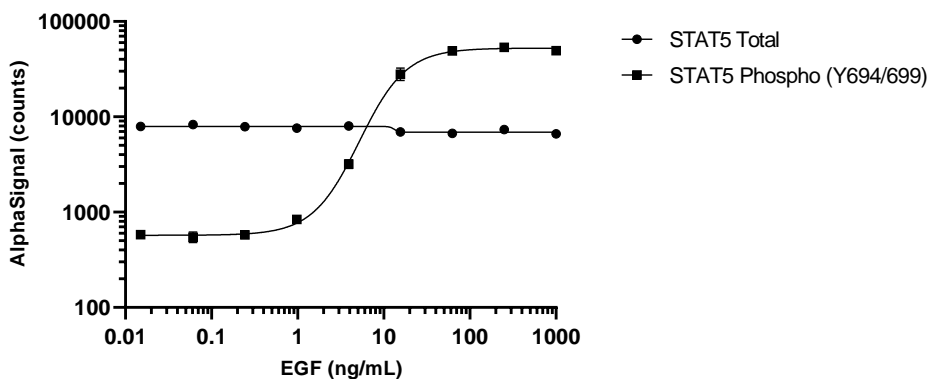
This assay kit contains antibodies which recognize the phospho-Tyr694 and Tyr699 epitopes, and a distal epitope, on STAT5A and STAT5B, respectively. The proteins detected by this kit correspond to UniProt ID P42229 and P51692. These antibodies recognize STAT5 of human origin. Other species should be 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 µg/mL EGF for 10 minutes and lysed in 8 mL of 1X *SureFire Ultra* Lysis Buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. A431 cells were plated at 40K cells per well in a 96 well plate and incubated overnight. Cells were serum starved for 2 hours prior to treatment with EGF at the indicated concentrations for 15 minutes. Cells were lysed in 1x Lysis Buffer and assayed separately for Total or Phospho (Tyr694/699) STAT5 using respective *SureFire Ultra* kits. Equivalent to approximately 4000 cells/datapoint.



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>