

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex p-STAT5 (Tyr694/695) + Total STAT5 Assay Kit

**Products: MPSU-PTST5-K500, MPSU-PTST5-K10K, MPSU-PTST5-K50K,
MPSU-PTST5-K-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Tyr694/695 epitope, and a distal epitope, on STAT5 for phospho STAT5 analysis, plus antibodies to measure the total levels of STAT5. The proteins detected by this kit correspond to GenBank Accessions NP_003143 and NP_036580. STAT5A is also known as MGF and STAT5; STAT5B is also known as STAT5.

These antibodies recognize STAT5 of human and mouse origin. They are predicted to react also with rat. Other species should be tested on a case-by-case basis.

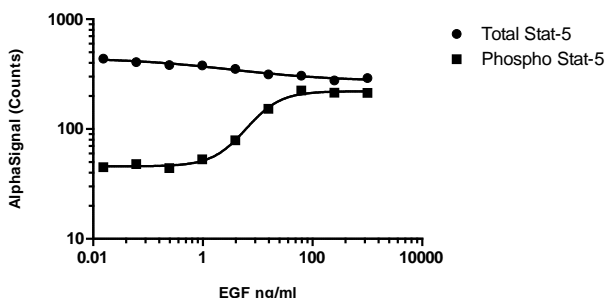
Control Lysate Information

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 8mL 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 15min, then lysed with 1X Lysis buffer and assayed for both phospho and total STAT5 as detailed in the MPSU manual.

EGF Dose Response in A431 cells - Stat-5 MPSU assay



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>