

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ STAT3 Total Assay Kit

**Products: ALSU-TST3-A500, ALSU-TST3-A10K, ALSU-TST3-A50K,
ALSU-TST3-A-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize two distinct invariant epitopes on STAT3 α and STAT3 β . The proteins detected by this kit correspond to GenBank Accessions NP_644805 and NP_998827.

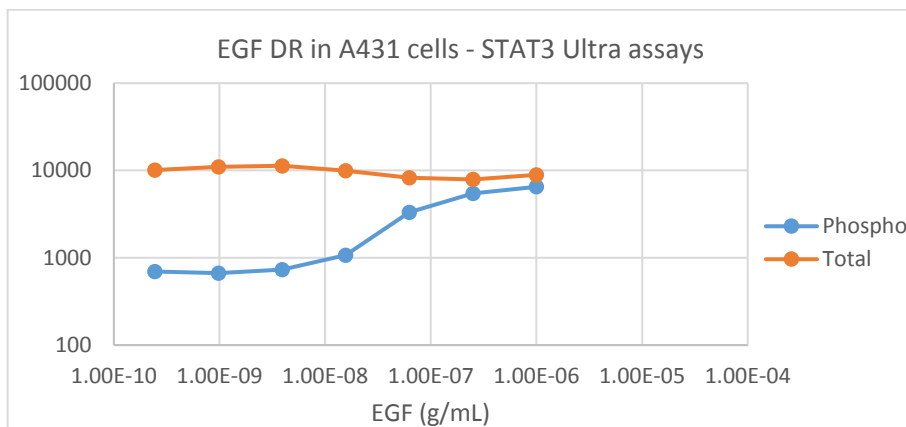
These antibodies recognize STAT3 of human, mouse and rat origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from A431 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 2 μ g/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. Lysates were assayed for Phospho and Total STAT3 in separate wells using the respective *SureFire Ultra* kits.



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>