

Technical Data Sheet

Caution: Research use only.

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

**Products: ALSU-PST5-B500, ALSU-PST5-B10K, ALSU-PST5-B50K,
ALSU-PST5-B-HV**

Kit-Specificity Information

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 GenBank Accessions NP_003143 and NP_036580. STAT5A is also known as MGF and STAT5; STAT5B is also known as STAT5.

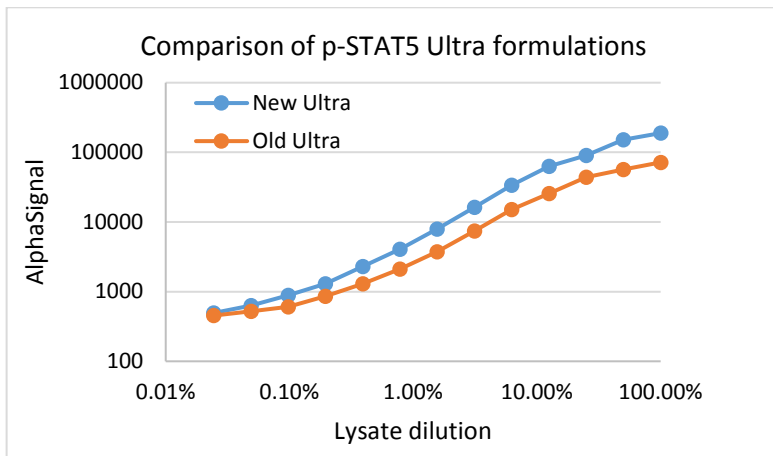
These antibodies recognize STAT5 of human origin. Other species should be tested on a case-by-case basis.

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 8mL of 1X *SureFire Ultra* Lysis buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. Data is shown for this new formulation kit and the previous old formulation kit.



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>
TGRSU196.0