

Alpha SureFire® Ultra™ Multiplex

Human p-STAT5 (Tyr694/699) + Total STAT5 Detection Kit

Product number: MPSU-PTST5-K500, MPSU-PTST5-K10K,
MPSU-PTST5-K50K, MPSU-PTST5-K-HV



Kit specificity:

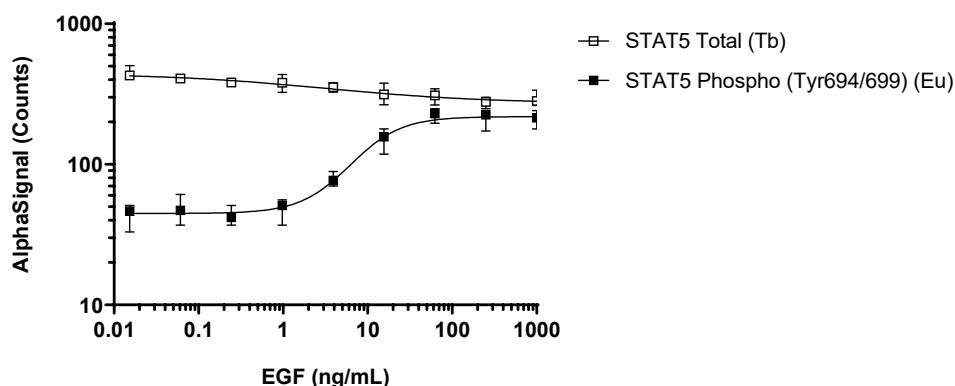
This assay kit contains antibodies which recognize the phospho-Tyr694/695 epitope and a distal epitope on STAT5, plus antibodies to measure the total levels of STAT5. The proteins detected by this kit correspond to UniProt IDs P42229 (STAT5A) and P51692 (STAT5B). STAT5 is also known as Signal Transducer and Activator of Transcription 5. 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 media, treated with 2 µg/mL rhEGF for 10 minutes and lysed with 8 mL of Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. A431 seeded at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with EGF at the indicated concentrations for 15 minutes. Cells were lysed with Lysis buffer and assayed for both Phospho (Tyr694/699) and Total STAT5 in the same wells of an assay plate using the SureFire Ultra Multiplex kit. Equivalent to approximately 4,000 cells/datapoint.



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF"