

AlphaLISA® SureFire® Ultra™ Multiplex

Human p-STAT6 (Tyr641) + Total STAT6 Detection Kit

Product number: MPSU-PTST6-K500, MPSU-PTST6-K10K,
MPSU-PTST6-K50K, MPSU-PTST6-K-HV



Kit specificity:

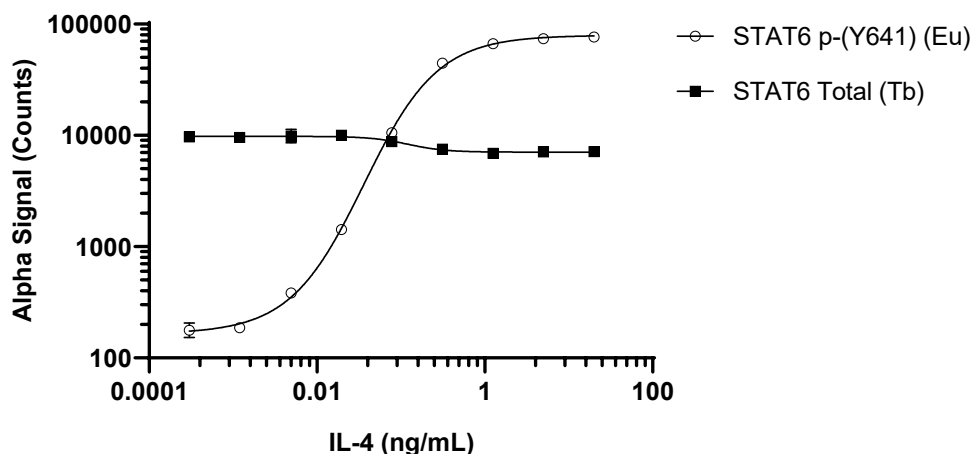
This assay kit contains antibodies which recognize the phospho-Tyr641 epitope and a distal epitope on STAT6 plus antibodies to detect Total STAT6. The proteins detected by this kit correspond to UniProt ID P42226. STAT6 is also known as IL-4 Stat. These antibodies recognize STAT6 of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

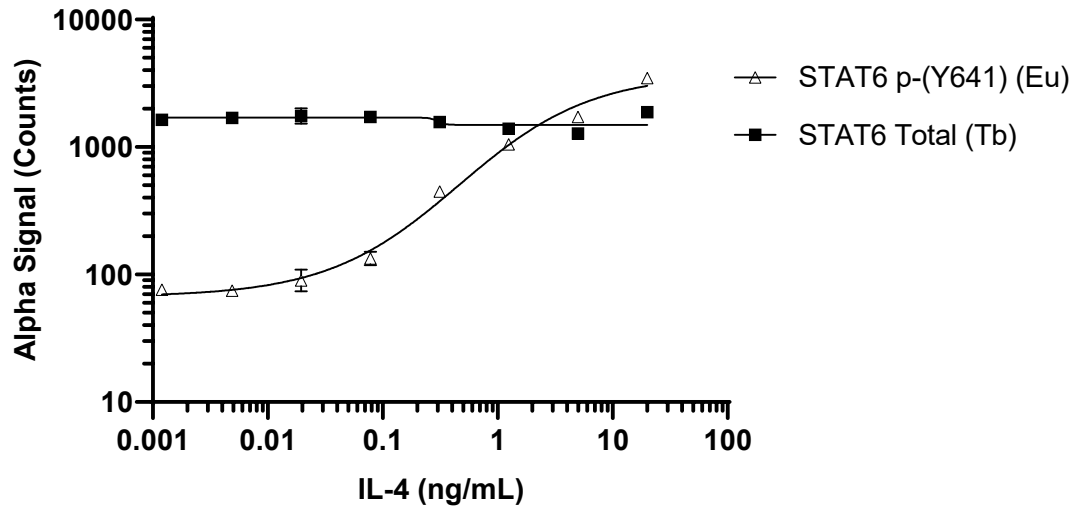
Positive Control Lysate: Prepared from HeLa cells, cultured to confluence in T175 flasks in 10% FBS containing medium. Cells were treated with 20 ng/mL recombinant human IL-4 for 20 minutes and lysed with 4 mL of Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. A431 cells were seeded at 50K cells/well in a 96 well plate and incubated overnight. Cells were treated with recombinant human IL-4 for 20 minutes. Cells were lysed with Lysis Buffer and assayed for Phospho (Tyr641) and Total STAT6 in the same wells of an assay plate using the SureFire Ultra Multiplex kit. Equivalent to approximately 2,500 cells/datapoint.



Data obtained with a 2-plate, 2-incubation protocol. THP-1 cells were harvested, then washed and resuspended in HBSS + 0.1 % BSA at 400K cells/well in a 96 well plate. Cells were treated with recombinant human IL-4 for 20 minutes. Cells were lysed with 5X Lysis Buffer and assayed for Phospho (Tyr641) and Total STAT6 in the same wells of an assay plate using the *SureFire Ultra* Multiplex kit. Equivalent to approximately 12,000 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"

