

AlphaLISA™ SureFire® Biotin Free

Human and Mouse STAT6 Total Detection Kit

Product number: ASBF-TST6-A500, ASBF-TST6-A10K,
ASBF-TST6-A50K, ASBF-TST6-A-HV



Kit specificity:

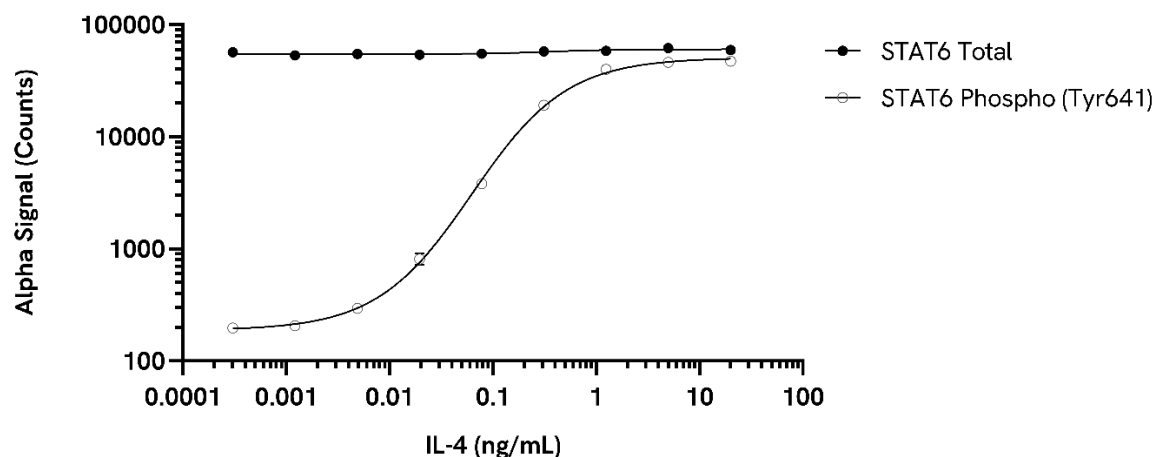
This assay kit contains antibodies which recognize distinct epitopes on signal transducer and activation of transcription 6 (STAT6). The protein detected by this kit corresponds to UniProt ID P42226. STAT6 is also known as IL-4 Stat. These antibodies recognize STAT6 of human and mouse origin. Other species 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, then treated with 20 ng/mL IL-4 for 20 minutes and lysed with 4 mL of Lysis Buffer.

Representative data:

Data obtained with a 2-plate, 2-incubation protocol. THP-1 cells were seeded at 400K cells/well in a 96 well plate in complete RPMI 1640 medium and treated with IL-4 at the indicated concentrations for 20 minutes. Cells were lysed with the addition of 5X Lysis Buffer and assayed separately for Phospho (Tyr641) and Total STAT6 using respective SureFire Biotin Free kits. Equivalent to approximately 16,000 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"