

AlphaLISA® SureFire® Ultra™

Human STAT4 Total Detection Kit

Product number: ALSU-TST4-A-500, ALSU-TST4-A-10K,
ALSU-TST4-A-50K, ALSU-TST4-A-HV



Kit specificity:

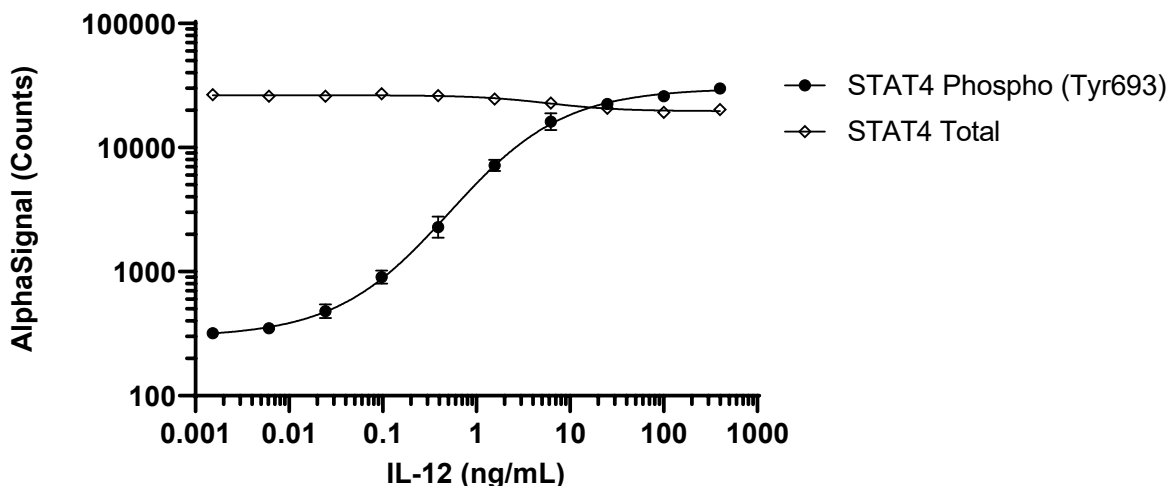
This assay kit contains antibodies which recognize distinct epitopes on signal transducer and activator of transcription 4 (STAT4). The protein detected by this kit corresponds to UniProt ID Q14765. These antibodies recognize STAT4 of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

Positive Control Lysate: Prepared from NK-92 cells, cultured to confluence in T175 flasks with MEM- α media containing 12.5% FBS, 12.5% Horse Serum, 0.2 mM inositol, 0.1 mM 2-mercaptoethanol, 0.02 mM folic acid and 200 U/mL recombinant IL-2. Cells were then deprived of IL-2 for 18 hours, treated with 200 ng/mL IL-12 for 30 minutes and lysed at 3.2×10^6 cells/mL with Lysis Buffer.

Representative data:

Data obtained with a 2-plate, 2-incubation protocol. NK-92 cells were seeded in a T175 flask at 200K/mL in media deprived of IL-2 and incubated overnight. Cells were washed in HBSS, seeded at 3×10^6 cells/mL in a 96 well plate, then treated with IL-12 at the indicated concentrations for 30 minutes. Cells were lysed with Lysis Buffer at 1.2×10^6 cells/mL and assayed separately for Phospho (Tyr693) and Total STAT4 using respective *SureFire Ultra* kits. Equivalent to 2,400 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"



Revvity, Inc.
940 Winter Street
Waltham, MA 02451 USA

(800) 762-4000
www.revvity.com

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2023, Revvity, Inc. All rights reserved.

VersionTGRSU278.3