

AlphaLISA® SureFire® Ultra™

# Human p-STAT4 (Tyr693) Detection Kit

**Product number:** ALSU-PST4-A500, ALSU-PST4-A10K,  
ALSU-PST4-A50K, ALSU-PST4-A-HV



**Kit specificity:**

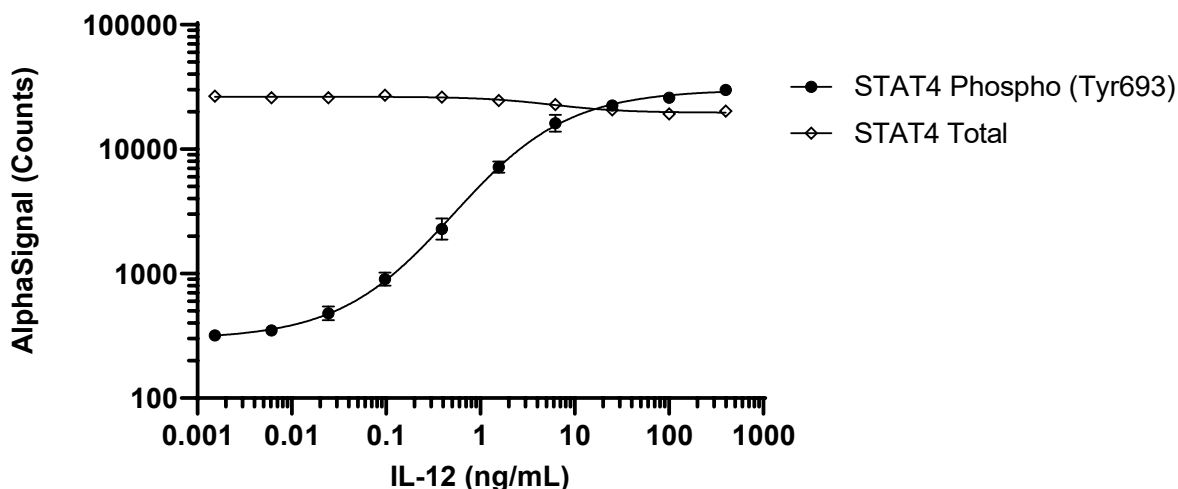
This assay kit contains antibodies which recognize the phospho-Tyr693 epitope and a distal epitope on signal transducer and activator of transcription 4 (STAT4). The protein detected by this kit corresponds to UniProt ID Q14765. These antibodies recognize STAT4 of human origin. Other species should be 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- $\alpha$  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  $3.2 \times 10^6$  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  $3 \times 10^6$  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  $1.2 \times 10^6$  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](http://www.revvity.com)

For a complete listing of our global offices, visit [www.revvity.com](http://www.revvity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.

VersionTGRSU195.3