

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ STAT4 Total Assay Kit

ALSU-TST4-A500, ALSU-TST4-A10K, ALSU-TST4-A50K, ALSU-TST4-A-HV

Kit-Specificity

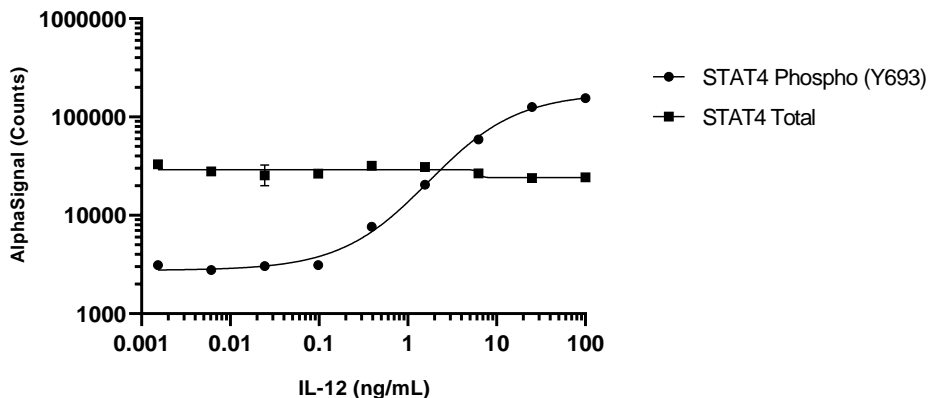
This assay kit contains antibodies which recognize two invariant epitopes on STAT4. The protein detected by this kit corresponds to UniProt ID Q14765. STAT4 is also known as Signal transducer and activator of transcription 4. These antibodies recognize STAT4 of human origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from NK92 cells, cultured to 0.2×10^6 cells/mL in T175 flasks in 12.5% FBS containing medium, IL-2 starved, then treated with 200 ng/mL IL-12 for 30 minutes and lysed at 3.2×10^6 cells/mL with 1x *SureFire Ultra* Lysis Buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. NK-92 cells were seeded in a flask at a density of 200K cells/mL in IL-12 free media and incubated overnight. Cells were then washed in HBSS and seeded at 8×10^6 cells/mL in a 96 well plate and then treated with IL-12 at the indicated concentrations for 30 minutes. Cells were lysed with 5x Lysis Buffer at a final density of 3.2×10^6 cells/mL and assayed for Total or Phospho (Tyr693) STAT4 using respective *SureFire Ultra* kits. Equivalent to 32 000 cells /datapoint.



For more information and detailed protocols, please refer to the generic AlphaLISA *SureFire Ultra* Assay Kits Manual, available from <http://www.perkinelmer.com/category/alpha-surefire-kits>