

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex

Terbium SureFire Ultra p-STAT4 (Tyr693) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

Products: TBSU-PST4-A500, TBSU-PST4-A10K, TBSU-PST4-A50K, TBSU-PST4-A-HV

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Tyr693 epitope, and a distal epitope, on STAT4. The protein detected by this kit corresponds to Entrez-Gene Id 6775. These antibodies recognize STAT4 of human origin. It is predicted to react also with mouse. Other species should be tested on a case-by-case basis.

The antibodies in this kit are identical to those in the corresponding ALSU-PST4-A kits, other than the CaptSure2 tag.

Control Lysate Information

Positive Control Lysate: Prepared from NK92 cells treated with IL-12 (50ng/mL) and lysed in 1X SureFire Ultra Lysis buffer (1x10⁶ cells/mL final lysate concentration).

Representative data

Control lysate samples measured in the terbium channel using an Envision plate reader.

