

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex

Terbium SureFire Ultra p-STAT1 (Ser727) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

**Products: TBSU-PST1-B500, TBSU-PST1-B10K, TBSU-PST1-B50K,
TBSU-PST1-B-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Ser727 epitope, and a distal epitope, on signal transducer and activator of transcription 1 (STAT1). The protein detected by this kit corresponds to GenBank Accession NP_009330. STAT1 is also known as ISGF-3, STAT1 and DKFZp686B04100. These antibodies recognize STAT1 of human and mouse origin. Other species should be tested on a case-by-case basis.

The antibodies in this kit are identical to those in the corresponding ALSU-PST1-B kits, other than the CaptSure2 tag.

Control Lysate Information

Positive Control Lysate: Prepared from Hela cells, cultured to confluence in T175 flasks in MEM with 10% FBS for 5 days, then treated with sorbitol (0.6M) for 45min, and lysed in 6mL of 1X SureFire Ultra Lysis buffer.

Representative data

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

