

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex

Terbium SureFire Ultra p-STAT1 (Tyr701) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

Products: TBSU-PST1-A500, TBSU-PST1-A10K, TBSU-PST1-A50K, TBSU-PST1-A-HV

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Tyr701 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-A 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 20min, and then stimulated with gamma-Interferon (200ng/mL) for 20 min, 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.

