

# Technical Data Sheet

Caution: Research use only.

## Alpha SureFire® Ultra™ Multiplex

### Terbium SureFire Ultra p-STAT5 (Tyr694/699) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

**Products: TBSU-PST5-B500, TBSU-PST5-B10K, TBSU-PST5-B50K,  
TBSU-PST5-B-HV**

#### **Kit-Specificity Information**

This assay kit contains antibodies which recognize the phospho-Tyr694 and Tyr699 epitopes, and a distal epitope, on STAT5A and STAT5B, respectively. The proteins detected by this kit correspond to GenBank Accessions NP\_003143 and NP\_036580. STAT5A is also known as MGF and STAT5; STAT5B is also known as STAT5. These antibodies recognize STAT5 of human 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-PST5-B kits, other than the CaptSure2 tag.

#### **Control Lysate Information**

Positive Control Lysate: Prepared from A431 cells, cultured to confluence in T175 flasks in 10% FBS containing medium for 3 days, then treated with EGF (2µg/mL) for 10min and lysed in 13mL of 1X SureFire Ultra Lysis buffer.

#### **Representative data**

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

