

# Technical Data Sheet

Caution: Research use only.

## Alpha SureFire® Ultra™ Multiplex

### Terbium SureFire Ultra p-STAT3 (Tyr705) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

**Products: TBSU-PST3-A500, TBSU-PST3-A10K, TBSU-PST3-A50K,  
TBSU-PST3-A-HV**

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

This assay kit contains antibodies which recognize the phospho-Tyr705-epitope, and a distal epitope on STAT3 $\alpha$  and STAT3 $\beta$ . The proteins detected by this kit correspond to GenBank Accessions NP\_644805 and NP\_998827. These antibodies recognize STAT3 of human, mouse and rat 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-PST3-A 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 (2 $\mu$ g/mL) for 10min and lysed in 10mL of 1X SureFire Ultra Lysis buffer.

#### **Representative data**

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

