

# Technical Data Sheet

Caution: Research use only.

## Alpha SureFire® Ultra™ Multiplex

### Terbium SureFire Ultra p-STAT6 (Tyr641) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

**Products: TBSU-PST6-A500, TBSU-PST6-A10K, TBSU-PST6-A50K, TBSU-PST6-A-HV**

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

This assay kit contains antibodies which recognize the phospho-Tyr641 epitope, and a distal epitope, on STAT6. These antibodies recognize STAT6 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-PST6-A kits, other than the CaptSure2 tag.

#### **Control Lysate Information**

Positive Control Lysate: Prepared from confluent HeLa cells in a T175 flask stimulated with 20 ng/mL IL-4 for 20 mins. Cells were lysed in 10 mL of 1X SureFire Ultra Lysis buffer.

#### **Representative data**

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

