

# Technical Data Sheet

Caution: Research use only.

## Alpha SureFire® Ultra™ Multiplex

### Terbium SureFire Ultra p-SYK (Tyr525/526) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

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

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

This assay kit contains antibodies which recognize the phospho-Tyr525/526 epitope, and distal invariant epitopes, on SYK. The protein detected by this kit corresponds to Entrez-Gene ID 6850. These antibodies recognize SYK of human origin. It is predicted to react also with mouse and rat. Other species should be tested on a case-by-case basis. The antibodies in this kit are identical to those in the corresponding ALSU-PSYK-A kits, other than the CaptSure2 tag.

#### **Control Lysate Information**

Positive Control Lysate: Prepared from U937 cells seeded at  $2 \times 10^6$  cells/mL and treated with 100 mM pervanadate for 30 min in HBSS + 0.5%BSA, and then lysed with 5X SureFire Ultra Lysis buffer (1:5) to give a final cell concentration of  $10^6$  cells/mL.

#### **Representative data**

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

