

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex

Terbium SureFire Ultra p-SLP-76 (Ser376) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

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

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Ser376 epitope, and a distal epitope, on SLP-76. The protein detected by this kit corresponds to GenBank Accession U20158.1. These antibodies recognize SLP-76 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-PSLP-A kits, other than the CaptSure2 tag.

Control Lysate Information

Jurkat cells were grown in RPMI with 10% FBS, spun down and reconstituted in HBSS at 4,000,000 cells per mL, and stimulated with 11 mM H₂O₂ + 50 μM calyculin for 10 mins. Cells were then lysed with SureFire Ultra 5X lysis buffer (1:5 in culture medium) to a final cellular density of 800,000 cells/mL.

Representative data

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

