

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex

Terbium SureFire Ultra Total Cofilin Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

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

Kit-Specificity Information

This assay kit contains antibodies which recognize two invariant epitopes on Cofilin. The protein detected by this kit corresponds to GenBank Accession NP_005498.1. Cofilin is also known as CFL and CFL1. These antibodies recognize Cofilin 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-TCOF-A kits, other than the CaptSure2 tag.

Control Lysate Information

Positive Control Lysate: Prepared from confluent HeLa cells grown in serum-containing MEM medium and lysed in 1X SureFire Ultra Lysis buffer.

Representative data

HeLa cells grown in 96 well plates were pretreated with AG1478 for 2 hours, then stimulated with EGF (1ng/mL) for 10 minutes, and cells lysed with 1X SureFire Ultra Lysis buffer. Lysates were then measured in multiplex mode for either p-ERK (Eu, ALSU kit) or Total Cofilin (Tb, TBSU kit) using an EnVision plate reader.

EGF stimulated Hela cells pretreated with AG1478

