

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex

Terbium SureFire Ultra p-IKK α (Ser176/180) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

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

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Ser176/180 epitope, and a distal epitope, on conserved helix-loop-helix ubiquitous kinase (IKK α). The protein detected by this kit corresponds to GenBank Accession NP_001269. IKK α is also known as IKK1, IKKA, IKBKA, TCF16, NFKB1A, IKK-alpha and CHUK. These antibodies recognize IKK α of human and 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-PIKKA-A kits, other than the CaptSure2 tag.

Control Lysate Information

Positive control: Prepared from recombinant p-IKK α protein at a concentration of 100 ng/mL in 1X Lysis Buffer. The controls are supplied lyophilized, and should be reconstituted in either dd H₂O or MilliQ® H₂O, and serially diluted as desired in 1X Lysis buffer. Once reconstituted, lysates should be stored frozen in single use aliquots.

Representative data

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

