

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex

Terbium SureFire Ultra p-NF-κB p65 (Ser536) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

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

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Ser536 epitope, and a distal epitope, on v-rel reticuloendotheliosis viral oncogene homolog A (avian) (NF-κB p65). The protein detected by this kit corresponds to GenBank Accession NP_068810. NF-κB p65 is also known as p65, NFKB3, MGC131774 and RELA. These antibodies recognize NF-κB p65 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-PNFKB-A kits, other than the CaptSure2 tag.

Control Lysate Information

Positive Control Lysate: Prepared from HeLa cells, cultured to confluence in T175 flasks in 10% FBS containing medium for 3 days, then treated with 50ng/mL TNFα and 20ng/mL Calyculin A for 10 minutes and lysed in 10mL of 1X SureFire Ultra Lysis buffer.

Representative data

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

