

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex p-NF-κB p65 (Ser536) + Total NF-κB p65 Assay Kit

**Products: MPSU-PTNFKB-K500, MPSU- PTNFKB-K10K, MPSU- PTNFKB-K50K,
MPSU- PTNFKB-K-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), plus antibodies to measure the total levels of 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.

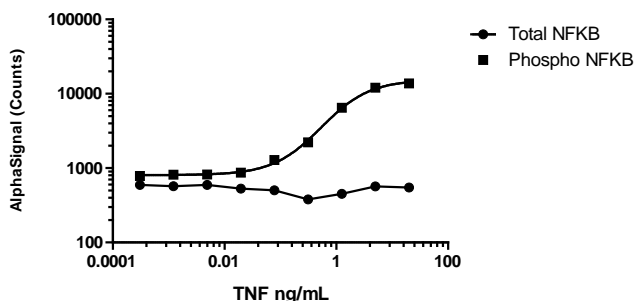
Control Lysate Information

Positive Control Lysate: Prepared from HeLa cells, cultured to confluence in T175 flasks in 10% FBS containing medium, 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

Data obtained with a 2-plate, 2-incubation protocol. Confluent HeLa cells were serum starved for 2 hours, then treated with TNFα at the indicated concentrations for 5min. Cells were then lysed with 1X Lysis buffer and assayed for both phospho and total NF-κB p65 as detailed in the MPSU manual.

TNF Dose Response in Hela cells - NF-κB MPSU



For more information and detailed protocols, please refer to the generic AlphaLISA *SureFire Ultra* Assay Kits Manual, available from <http://www.perkinelmer.com/category/alpha-surefire-kits>