

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ NFκB p65 Total Assay Kit

**Products: ALSU-TNFKB-A500, ALSU-TNFKB-A10K, ALSU-TNFKB-A50K,
ALSU-TNFKB-A-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize invariant epitopes on 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 NFκB p65 of human, rat and mouse origin. Other species should be tested on a case-by-case basis.

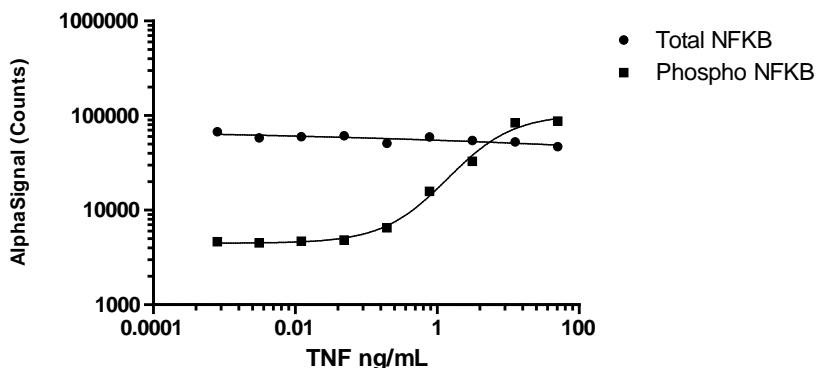
Control Lysate Information

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

Representative data

Data obtained with a 2-plate, 2-incubation protocol. MCF-7 cells were plated at 100K cells per mL, incubated overnight, starved for 2 hours, then stimulated with TNFα for 5 minutes at the indicated concentrations. Cells were lysed with 100μL of 1X Lysis buffer and assayed.

TNF Dose Response in Hela cells



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>.