

# Technical Data Sheet

## AlphaLISA® SureFire® Ultra™ NF-κB p65 Total Assay Kit

ALSU-TNFKB-A500, ALSU-TNFKB-A10K, ALSU-TNFKB-A50K, ALSU-TNFKB-A-HV

### Kit-Specificity

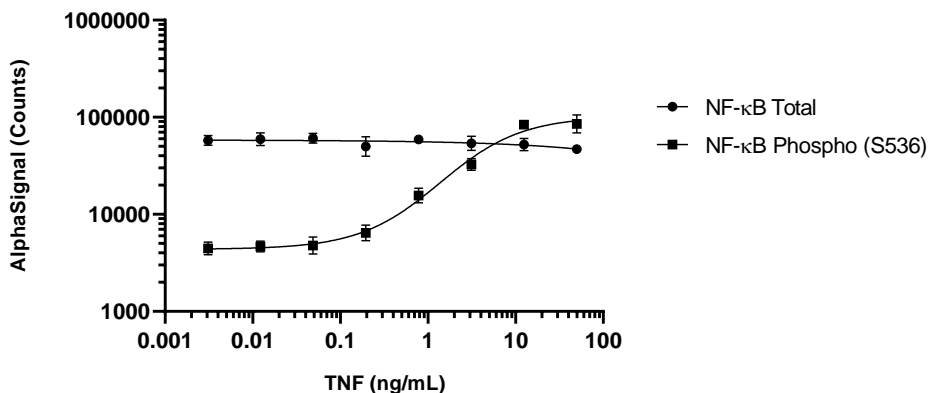
This assay kit contains antibodies which recognize invariant epitopes on NF-κB p65. The protein detected by this kit corresponds to UniProt ID Q04206. NF-κB p65 is also known as Transcription factor p65. These antibodies recognize NF-κB p65 of human, rat and mouse origin. Other species should be tested 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 50 ng/mL TNFα plus 20 ng/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. HeLa cells were plated at 40K cells per well in a 96 well plate and incubated overnight. Cells were starved for 2 hours and then stimulated with TNFα for 5 minutes at the indicated concentrations. Cells were lysed with 1X Lysis Buffer and assayed for Total or Phospho (Ser536) NF-κB, using respective *SureFire Ultra* assay kits. Equivalent to approximately 4000 cells/datapoint.



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>.