

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ p-NF-κB p65 (Ser536) Assay Kit

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

Kit-Specificity Information

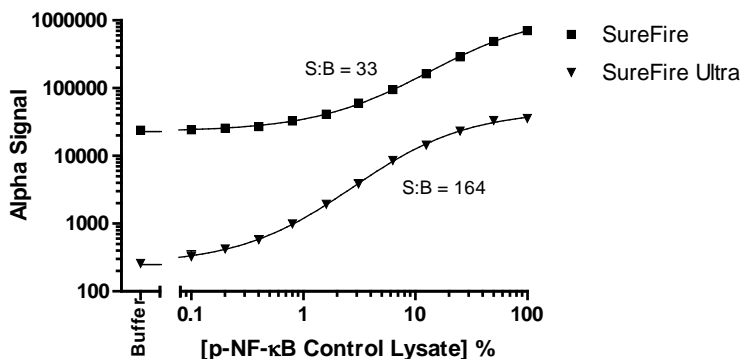
This assay kit contains antibodies which recognize the phospho-Ser536 epitope, and a distal epitope, on v-rel reticuloendotheliosis viral oncogene homolog A (avian) (NF-κB p65). The protein detected by this kit corresponds to GenBank Accession NP_068810. NF-κB p65 is also known as p65, NFKB3, MGC131774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

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. Data shows a comparison of AlphaScreen SureFire and AlphaLISA SureFire Ultra assay performance. Control lysate was diluted in the Lysis buffer specific for the assay type.



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>