

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ p-TBK1 (Ser172) Assay Kit

ALSU-PTBK1-A500, ALSU-PTBK1-A10K, ALSU-PTBK1-A50K, ALSU-PTBK1-A-HV

Kit-Specificity

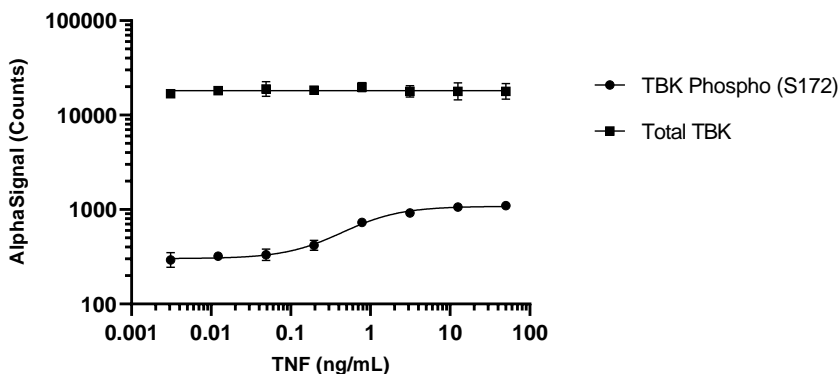
This assay kit contains antibodies that recognize the phospho Ser172 epitope and an invariant epitope on TBK1 (TANK-binding kinase 1). The protein detected by this kit corresponds to UniProt ID Q9UHD2. TBK1 is also known as NAK. These antibodies recognize TBK1 of human, mouse and rat origin. Other species 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 50 ng/mL TNF α and 20 ng/mL Calyculin A for 15 minutes, and lysed in 5 mL 1X *SureFire Ultra* Lysis Buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. HeLa cells plated in a 96 well plate overnight at 200K cells/mL were then treated with the indicated concentrations of TNF α for 15 minutes, lysed with 50 μ L of 1X Lysis Buffer. Samples were assayed separately for Total and Phospho (Ser172) TBK1, using respective *SureFire Ultra* kits. Equivalent to approximately 8000 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>.