

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ TBK1 Total Assay Kit

**Products: ALSU-TTBK1-A500, ALSU-TTBK1-A10K, ALSU-TTBK1-A50K,
ALSU-TTBK1-A-HV**

Kit-Specificity Information

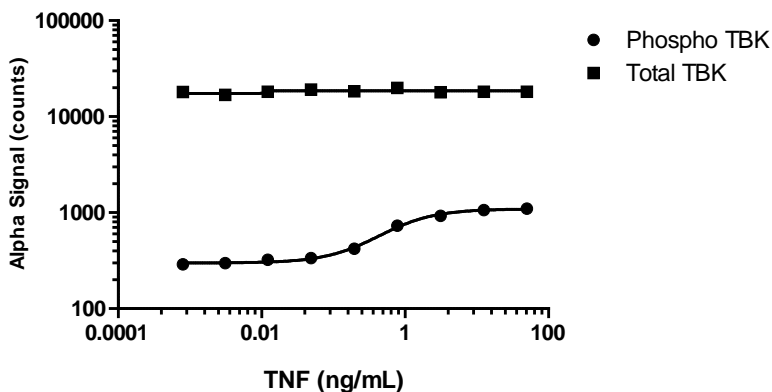
This assay kit contains antibodies that recognize invariant epitopes on TBK1 (TANK-binding kinase 1). The protein detected by this kit corresponds to Entrez-Gene ID 29110. TBK1 is also known as NAK. These antibodies recognize TBK1 of human, mouse and rat origin. Other species should be tested on a case-by-case 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 50ng/mL TNF α and 20ng/mL Calyculin A for 15 minutes, and lysed in 5ml 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.



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>.