

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ STING Total Assay Kit

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

Kit-Specificity

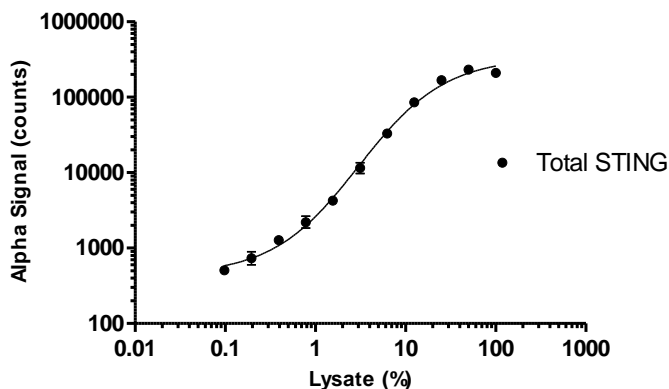
This assay kit contains antibodies which recognize invariant epitopes on Stimulator of interferon genes protein (STING). The protein detected by this kit corresponds to UniProt ID Q86WV6. STING is also known as ERIS, hMITA and transmembrane protein 173. These antibodies recognize STING of human origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from THP-1 cells, cultured at 500,000 cells/mL in T175 flasks in RPMI medium containing 10% FBS. Cells were harvested, washed with HBSS and lysed with 1x *SureFire Ultra* Lysis Buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. Control lysate (THP-1) was diluted with 1X Lysis Buffer and assayed for Total STING. The 100% value corresponds to approximately 8,000 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>.