

AlphaLISA® SureFire® Ultra™

Human STING Total Detection Kit

Product number: ALSU-TSTNG-A500, ALSU-TSTNG-A10K,
ALSU-TSTNG-A50K, ALSU-TSTNG-A-HV



Kit specificity:

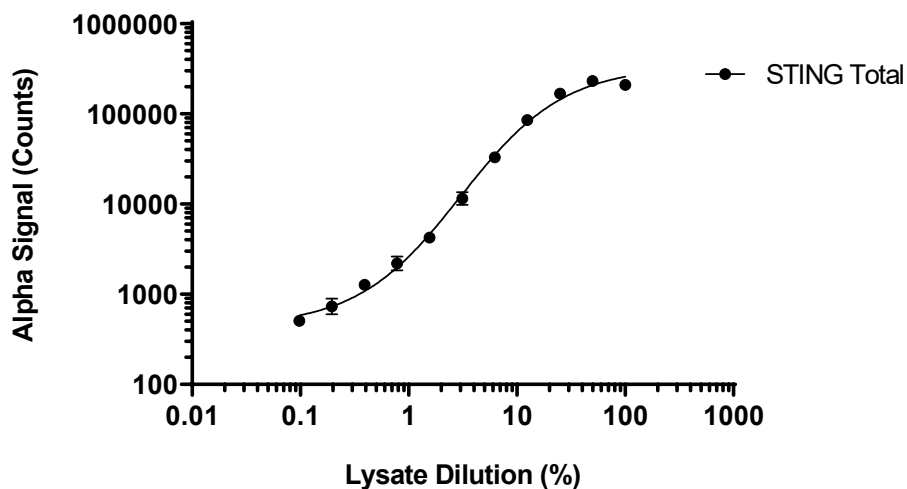
This assay kit contains antibodies which recognize distinct epitopes on Stimulator of interferon genes protein (STING). The protein detected by this kit corresponds to UniProt ID Q86WV6. STING is also known as ERIS and hMITA. These antibodies recognize STING of human origin. Other species should be tested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

Control lysate information:

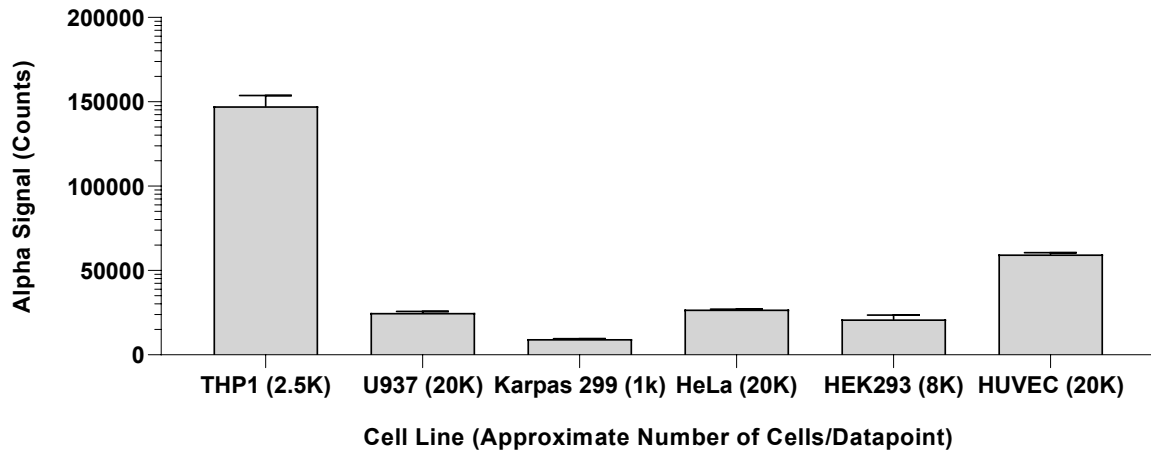
Positive Control Lysate: Prepared from THP-1 cells cultured in 10% FBS containing medium. Cells were washed in HBSS + 0.1% BSA, then spun down at 300 RCF for 5 minutes and lysed in Lysis Buffer at 1×10^6 cells/mL.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. Positive control lysate (THP-1) was serially diluted with Lysis Buffer and assayed for Total STING using the *SureFire Ultra* kit. At 100% lysate dilution, equivalent to approximately 8,000 cells/datapoint.



Data obtained from measurement of STING Total in various cell types lysed with Lysis Buffer. Approximate number of cell/datapoint is indicated for the various cell lines.



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF"

