

AlphaLISA® SureFire® Ultra™

Human TDP43 Total Detection Kit

Product number: ALSU-TTDP43-A500, ALSU TTDP43-A10K,
 ALSU-TTDP43-A50K, ALSU-TTDP43-A-HV



Kit specificity:

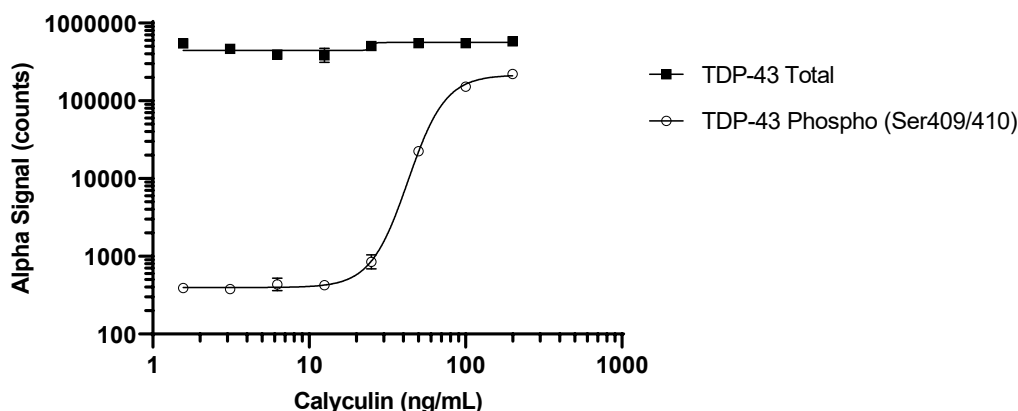
This assay kit contains antibodies which recognize distinct epitopes on TDP-43. The protein detected by this kit corresponds to UniProt ID Q13148. TDP-43 is also known as TAR DNA binding protein 43. These antibodies recognize TDP-43 of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

Positive Control Lysate: Prepared from U937 cells grown in 10% FBS containing medium. Cells were harvested, adjusted to 2×10^6 cells/mL and treated with 100 ng/mL Calyculin for 4 hours. Cells were spun down at 300 RCF for 5 minutes, washed with HBSS + 0.1% BSA and lysed with Lysis Buffer at 1×10^6 cells/mL.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. U937 cells were washed and resuspended in HBSS + 0.1% BSA and seeded at 200K cells/well in a 96 well plate. Cells were treated with Calyculin at the indicated concentrations for 4 hours. Cells were lysed with 5X Lysis Buffer and assayed separately for Phospho (Ser409/410) and Total TDP-43 using respective *SureFire Ultra* kits. Equivalent to approximately 8,000 cells/datapoint.



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"