

AlphaLISA® SureFire® Ultra™

Human PINK1 Total Detection Kit

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



Kit specificity:

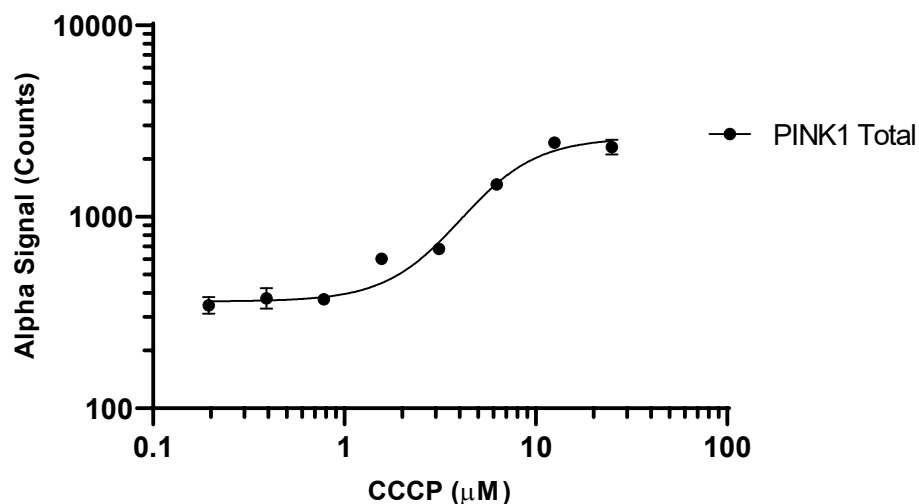
This assay kit contains antibodies which recognize distinct epitopes on PINK1. The protein detected by this kit corresponds to UniProt ID Q9BXM7. PINK1 is also known as PTEN induced putative kinase 1. These antibodies recognize PINK1 of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

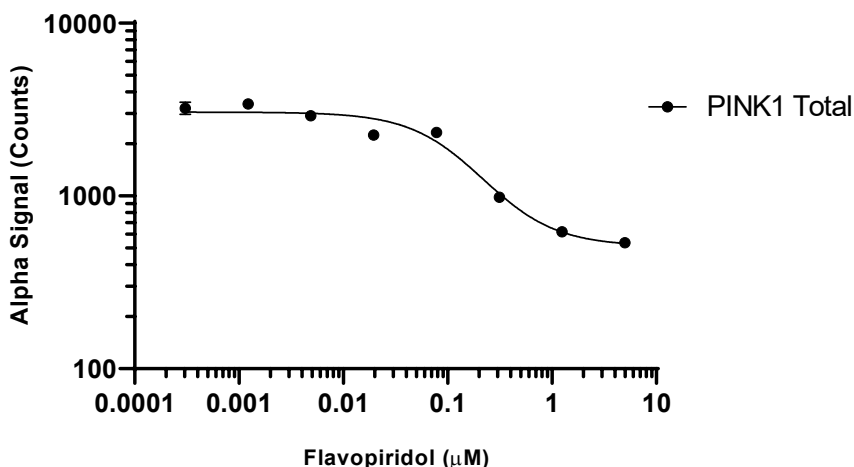
Positive Control Lysate: Prepared from PC-3 cells, cultured to confluence in T175 flasks in 10% FBS containing media, then treated with 25 μM CCCP for 18 hours and lysed with 4 mL of Lysis Buffer B.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. SHSY-5Y cells were plated at 10K cells/well in a 96 well plate and incubated for 48 hours. Cells were treated with CCCP at the indicated concentrations for 18 hours. Cells were lysed with Lysis Buffer B and assayed for Total PINK1 using the *SureFire Ultra* kit. Equivalent to approximately 4,000 cells/datapoint.



Data obtained with a 2-plate, 2-incubation protocol. HeLa cells were plated at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with Flavopiridol at the indicated concentrations in the presence of 12.5 μM CCCP for 16 hours. Cells were lysed with Lysis Buffer B and assayed for Total PINK1 using the *SureFire* Ultra kit. 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”

