

AlphaLISA® SureFire® Ultra™

Human AIOLOS Total Detection Kit

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



Kit specificity:

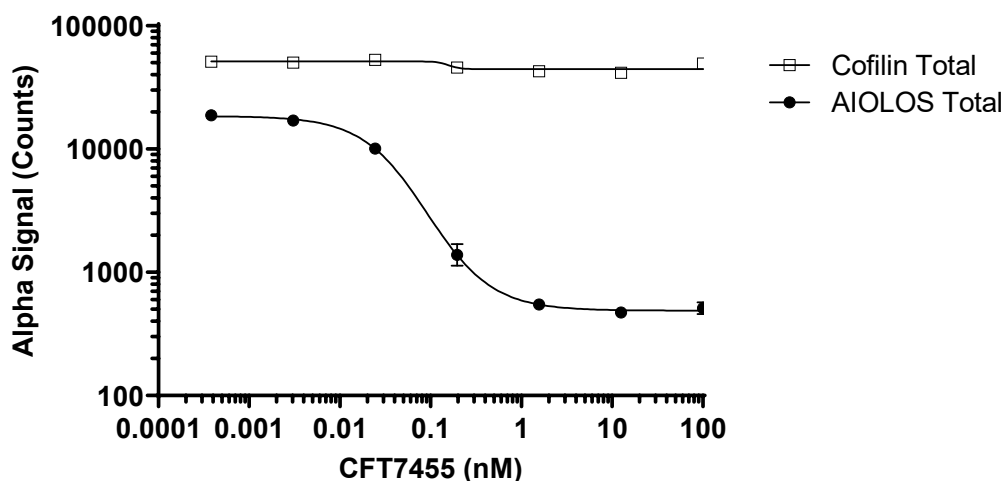
This assay kit contains antibodies which recognize distinct epitopes on AIOLOS. The protein detected by this kit corresponds to UniProt ID Q9UKT9. AIOLOS is also known as Zinc finger protein Aiolos and Ikaros family zinc finger protein 3. These antibodies recognize AIOLOS of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

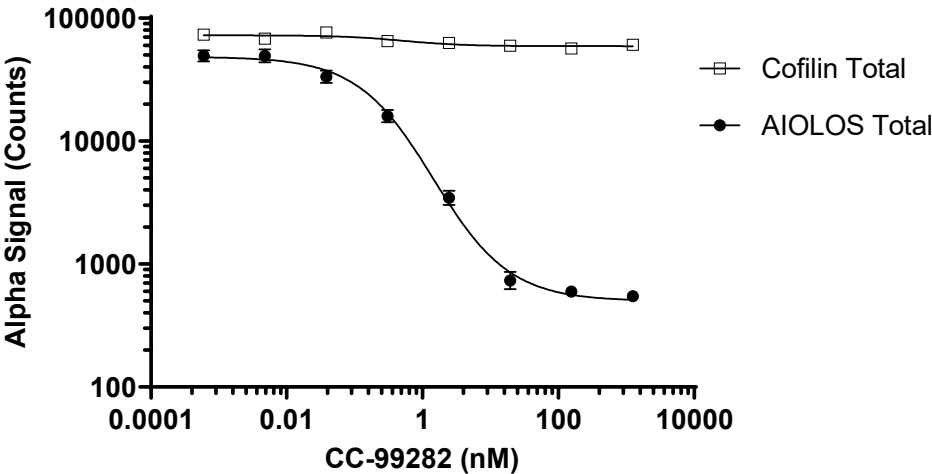
Positive Control Lysate: Prepared from Ramos cells, cultured to approximately 1×10^6 cells/mL in T175 flasks in 10% FBS containing medium and lysed at 5×10^6 cells/mL with Lysis Buffer.

Representative data

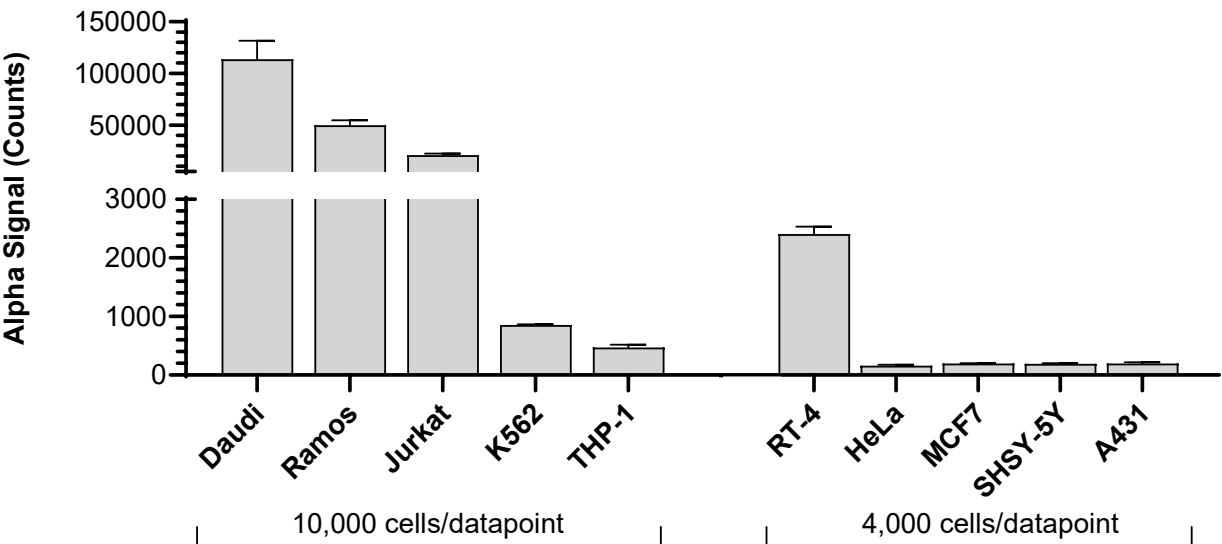
Data obtained with a 2-plate, 2-incubation protocol. Ramos cells were seeded at 100K cells/well in a round bottom 96 well plate. Cells were treated with CFT7455 at the indicated concentrations for 24 hours. Cells were lysed with Lysis Buffer and assayed separately for Total AIOLOS and Cofilin using respective SureFire Ultra kits. Equivalent to approximately 2,500 cells/datapoint (500 cells/datapoint for Cofilin).



Data obtained with a 2-plate, 2-incubation protocol. Ramos cells were seeded at 100K cells/well in a round bottom 96 well plate. Cells were treated with CC-99282 at the indicated concentrations for 24 hours. Cells were lysed with Lysis Buffer and assayed separately for Total AIOLOS and Cofilin using respective *SureFire Ultra* kits. Equivalent to approximately 2,500 cells/datapoint (500 cells/datapoint for Cofilin).



Data obtained from measurement of AIOLOS Total in various cell types lysed with Lysis Buffer. Approximate number of cells/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"

