

AlphaLISA® SureFire® Ultra™

Human and Mouse IKAROS Total Detection Kit

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



Kit specificity:

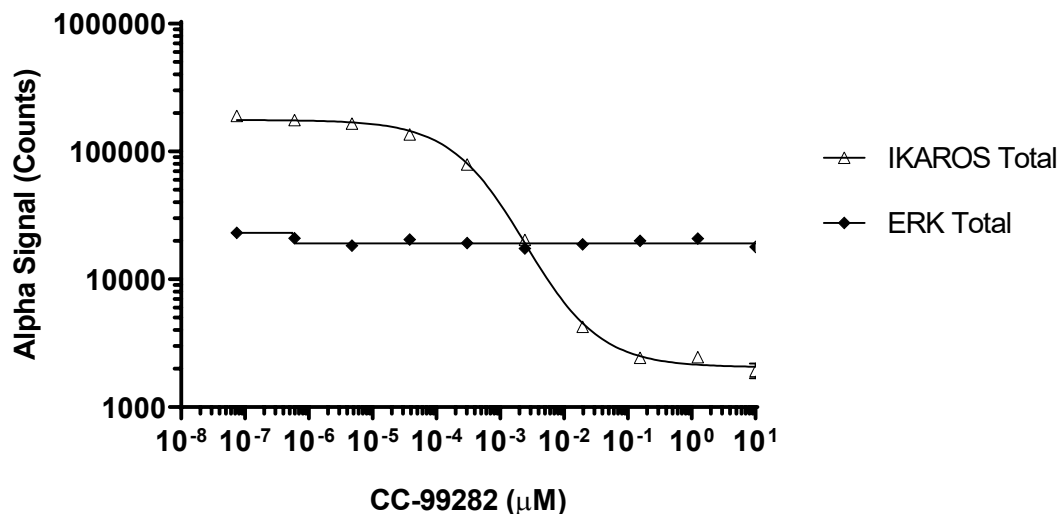
This assay kit contains antibodies which recognize distinct epitopes on IKAROS. The protein detected by this kit corresponds to UniProt ID Q13422. IKAROS is also known as Ikaros family zinc finger protein 1 (IKZF1), DNA-binding protein Ikaros, and Lymphoid transcription factor LyF-1. These antibodies recognize IKAROS of human and mouse origin. Other species should be tested on a case-by-case basis.

Control lysate information:

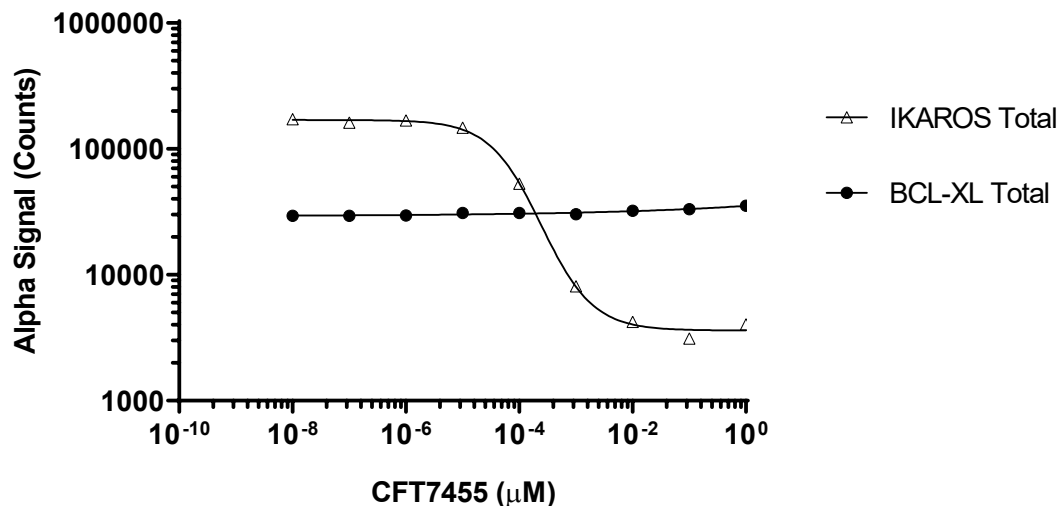
Positive Control Lysate: Prepared from Ramos cells, seeded at 500K/mL and grown overnight in T175 flasks in medium containing 10% FBS. Cells were washed in HBSS + 0.1% BSA and lysed in Lysis Buffer at a final cell concentration of 4×10^6 cells/mL.

Representative data

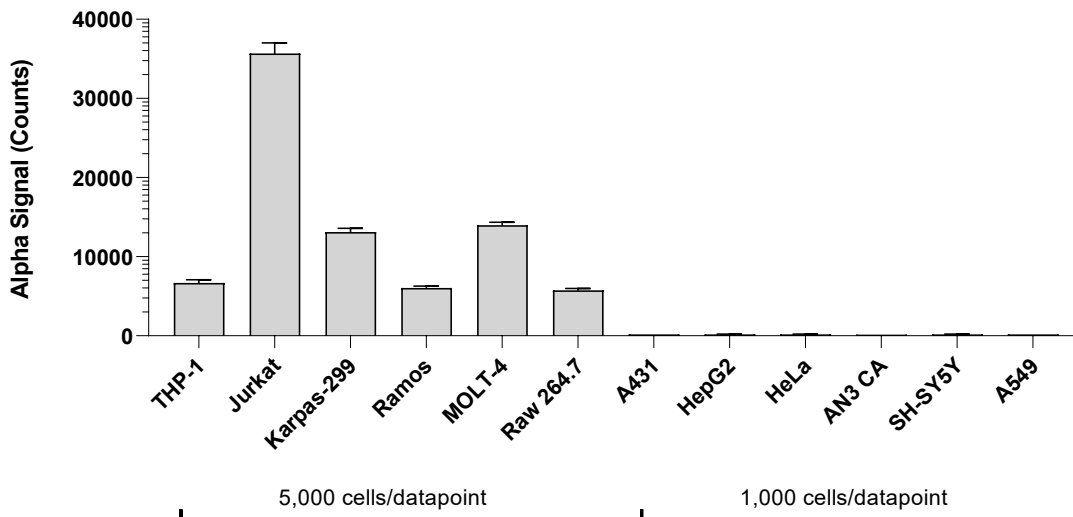
Data obtained with a 2-plate, 2-incubation protocol. Ramos cells were seeded at 200K/well in a 96 well plate and treated with CC-99282 at the indicated concentrations for 18 hours. Cells were washed with HBSS, lysed with Lysis Buffer and assayed separately for Total IKAROS and Total ERK using respective *SureFire Ultra* kits. Equivalent to approximately 10,000 cells/datapoint.



Data obtained with a 2-plate, 2-incubation protocol. MOLT-4 cells were seeded at 200K/well in a 96 well plate and treated with CFT7455 IKAROS PROTAC at the indicated concentrations for 18 hours. Cells were washed with HBSS, lysed with Lysis Buffer and assayed separately for Total IKAROS and Total BCL-XL using respective *SureFire Ultra* kits. Equivalent to approximately 10,000 cells/datapoint.



Data obtained for IKAROS Total from screening various lysates prepared from cells lysed in 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"

