

AlphaLISA® SureFire® Ultra™

Human and Mouse BCL-XL Total Detection Kit

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



Kit specificity:

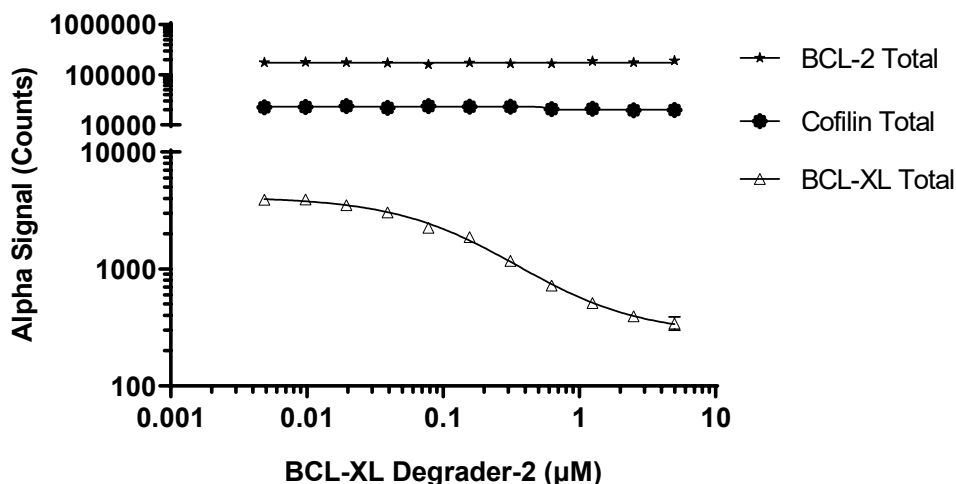
This assay kit contains antibodies which recognize distinct epitopes on BCL-XL. The protein detected by this kit corresponds to UniProt ID Q07817. BCL-XL is also known as Bcl-2-like protein 1, Bcl2-L-1 and Apoptosis regulator Bcl-X. These antibodies recognize BCL-XL of human and mouse origin. Other species should be tested on a case-by-case basis.

Control lysate information:

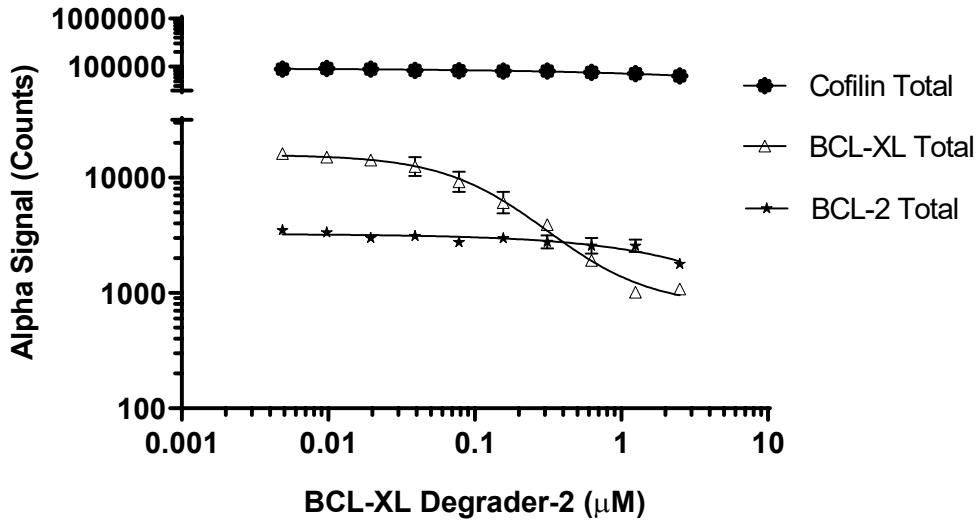
Positive Control Lysate: Prepared from A549 cells, cultured to confluence in T175 flasks in 10% FBS containing medium and lysed with 8mL of Lysis Buffer.

Representative data

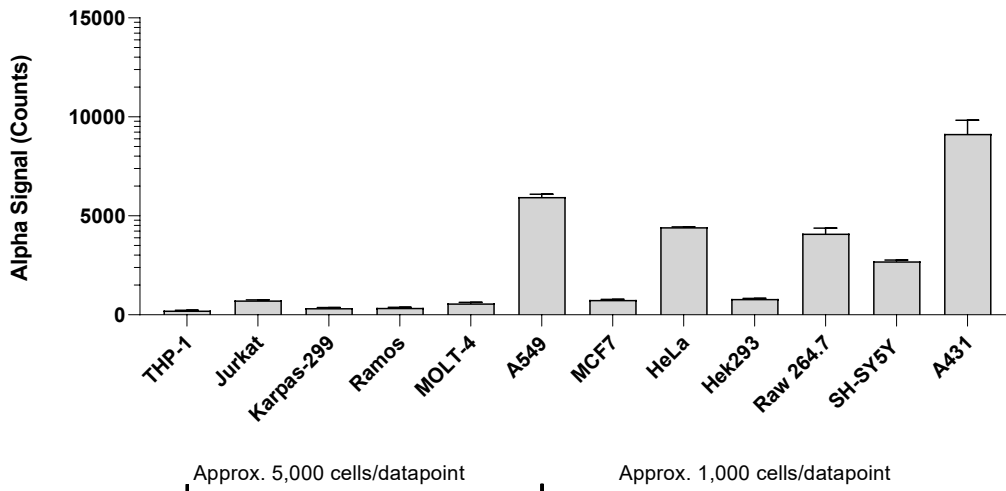
Data obtained with a 2-plate, 2-incubation protocol. SH-SY5Y cells were plated at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with BCL-XL Degradator 2 at the indicated concentrations for 22 hours. Cells were lysed with Lysis Buffer and assayed separately for BCL-2 Total, Cofilin Total and BCL-XL Total using respective SureFire Ultra kits. Equivalent to approximately 4,000 cells/datapoint or 80 cells/datapoint for Cofilin Total.



Data obtained with a 2-plate, 2-incubation protocol. Ramos cells were plated at 400K cells/well in a 96 well plate. Cells were treated with BCL-XL Degradator-2 at the indicated concentrations for 22 hours. Cells were washed with HBSS and lysed with Lysis Buffer and assayed separately for BCL-2 Total, Cofilin Total and BCL-XL Total using respective *SureFire Ultra* kits. Equivalent to approximately 20,000 cells/datapoint or 400 cells/datapoint for Cofilin Total.



Data obtained from measurement of BCL-XL 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"

