

AlphaLISA® SureFire® Ultra™

Human p-BCL-2 (Ser70) Detection Kit

Product number: ALSU-PBCL2-A500, ALSU-PBCL2-A10K,
ALSU-PBCL2-A50K, ALSU-PBCL2-A-HV



Kit specificity:

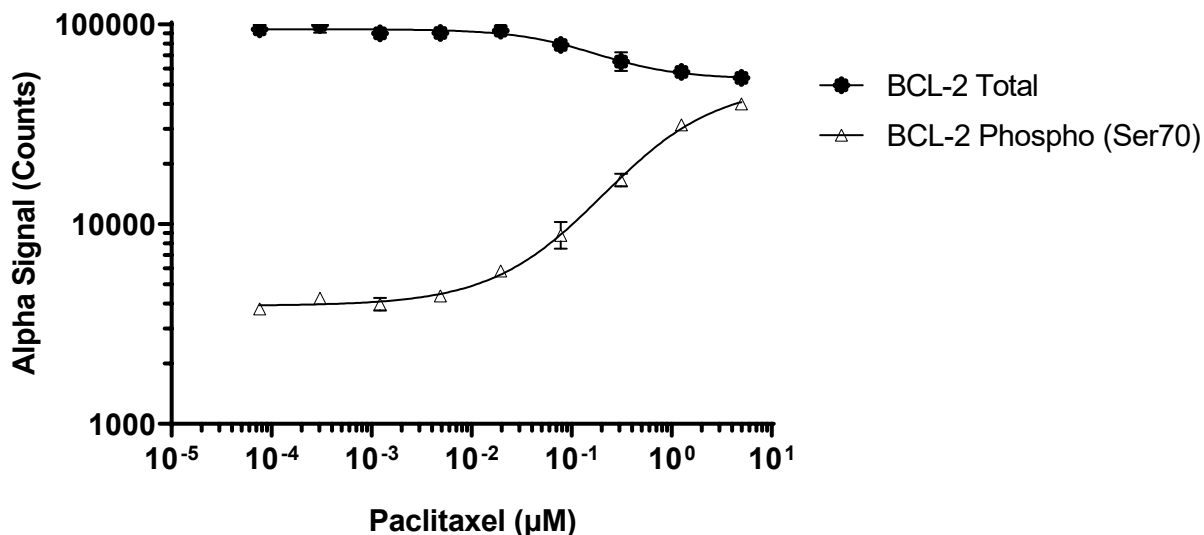
This assay kit contains antibodies which recognize the phospho-Ser70 epitope and a distal epitope on BCL-2. The protein detected by this kit corresponds to UniProt ID P10415. BCL-2 is also known as Apoptosis regulator Bcl-2. These antibodies recognize BCL-2 of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

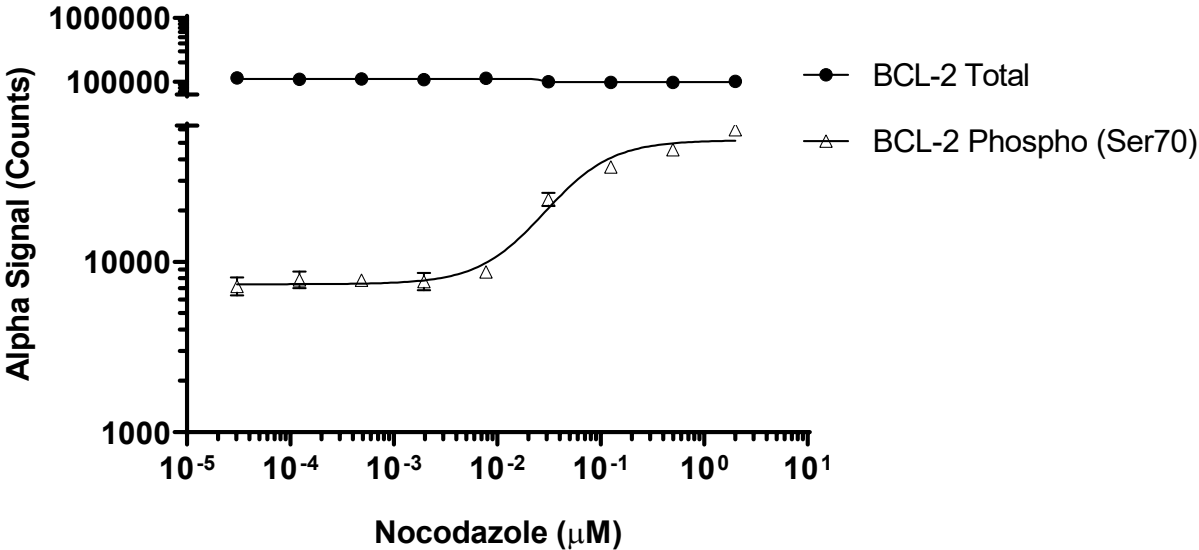
Positive Control Lysate: Prepared from SH-SY5Y cells, cultured to 90% confluence in T175 flasks in 10% FBS containing media, then treated with 1µM Paclitaxel for 24 hours and lysed with 5 mL Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. SH-SY5Y cells were plated at 20K cells/well in a 96 well plate and incubated overnight. Cells were treated with Paclitaxel at the indicated concentrations for 24 hours. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Ser70) and Total BCL-2 using respective *SureFire Ultra* kits. Equivalent to approximately 2,000 cells/datapoint.



Data obtained with a 2-plate, 2-incubation protocol. Jurkat cells were seeded at 200K cells/mL in a 96 well plate. Cells were treated with Nocodazole at the indicated concentrations for 18 hours. Cells were washed in HBSS and lysed in Lysis Buffer and assayed for Phospho BCL-2 (Ser70) and BCL-2 Total using respective *SureFire Ultra* kits. Equivalent to approximately 20,000 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"

