

AlphaLISA® SureFire® Ultra™

Human BRD4 Total Detection Kit

Product number: ALSU-TBRD4-A500, ALSU-TBRD4-A10K,
ALSU-TBRD4-A50K, ALSU-TBRD4-A-HV



Kit specificity:

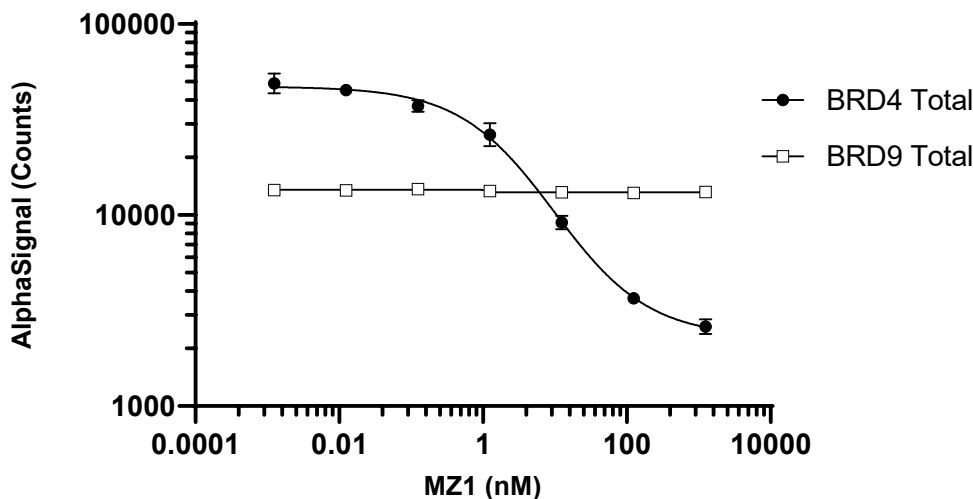
This assay kit contains antibodies which recognize distinct epitopes on BRD4. The protein detected by this kit corresponds to UniProt ID O60885. BRD4 is also known as Bromodomain-containing protein 4 and Protein HUNK1. These antibodies recognize BRD4 of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

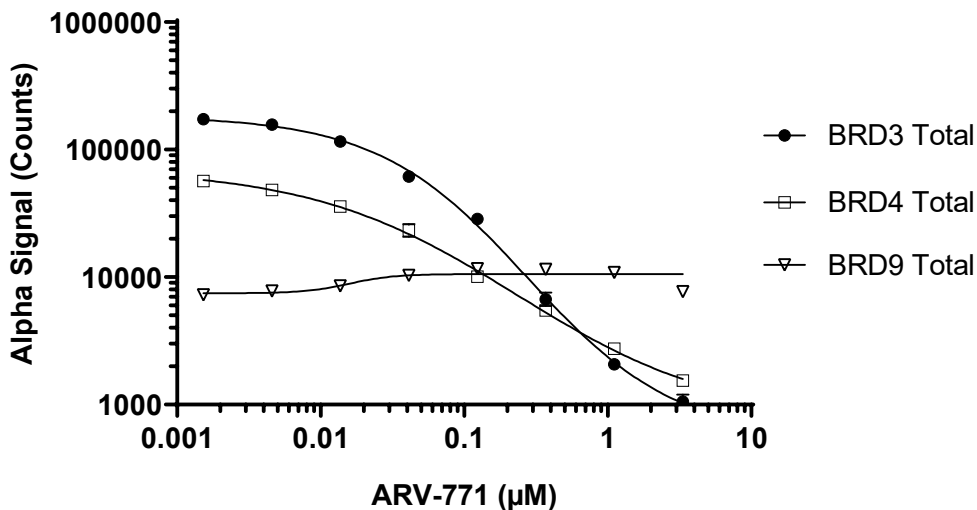
Positive Control Lysate: Prepared from RT4 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, lysed with 20mL of Lysis Buffer.

Representative data:

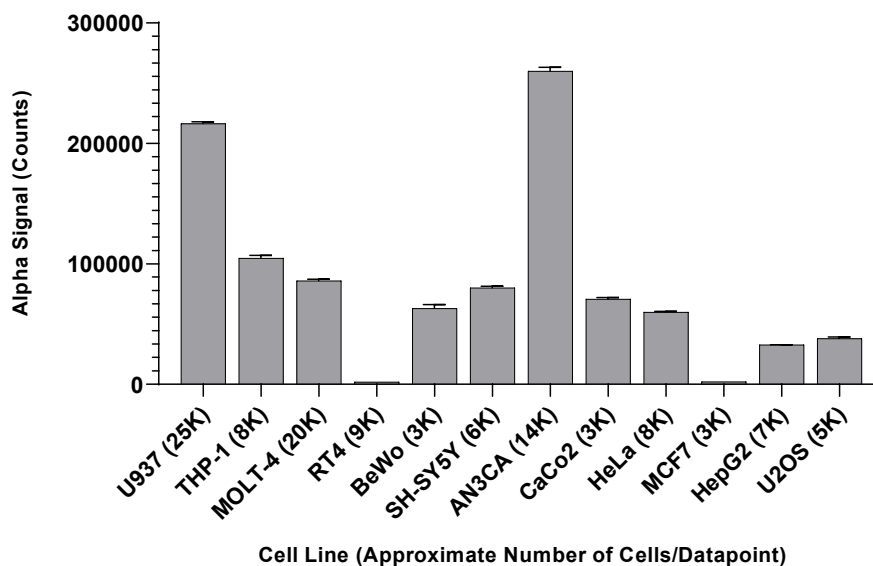
Data obtained with a 2-plate, 2-incubation protocol. THP-1 cells were harvested and seeded at 500K cells/well in a 96 well plate in HBSS containing 0.1% BSA. Cells were treated with MZ1 PROTAC at the indicated concentrations for 4 hours. Cells were lysed with the addition of 5X Lysis Buffer and lysates were assayed separately for BRD4 Total and BRD9 Total using respective Surefire Ultra kits. Equivalent to approximately 20,000 cells/datapoint.



Data obtained with a 2-plate, 2-incubation protocol. HEK293 cells were seeded at 20K cells/well in a 96 well plate and incubated overnight. Cells were treated with ARV-771 PROTAC at the indicated concentrations for 18 hours. Cells were lysed in Lysis Buffer and lysates were assayed separately for BRD4 Total, BRD3 Total and BRD9 Total using respective *Surefire Ultra* kits. Equivalent to approximately 2,000 cells/datapoint.



Data obtained with a 2-plate, 2-incubation protocol. Various cell types were lysed with Lysis Buffer and assayed for BRD4 Total using *SureFire Ultra* kit. 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"

