

AlphaLISA® SureFire® Ultra™

## Human WDR5 Total Detection Kit

**Product number:** ALSU-TWDR-A500, ALSU-TWDR-A10K,  
ALSU-TWDR-A50K, ALSU-TWDR-A-HV



### Kit specificity:

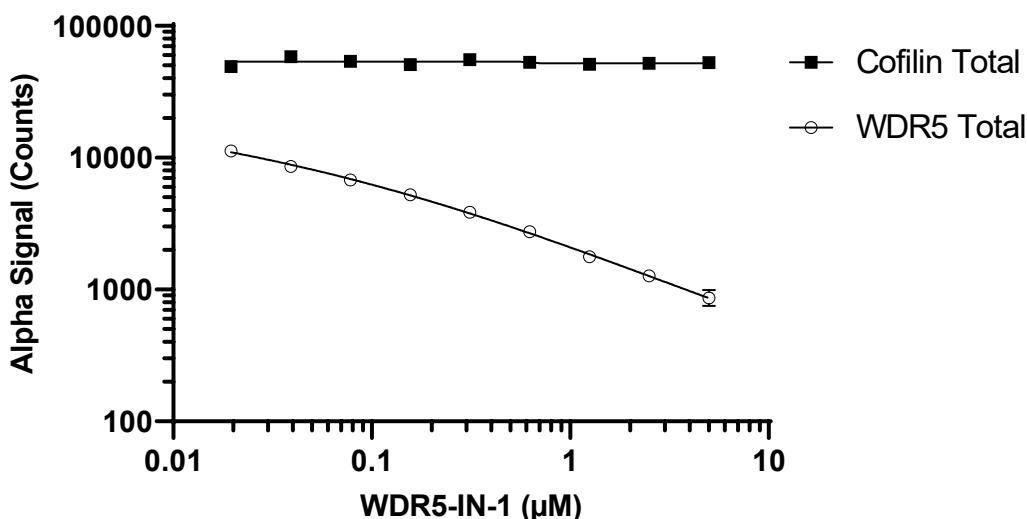
This assay kit contains antibodies which recognize distinct epitopes on WDR5. The protein detected by this kit corresponds to UniProt ID P61964. WDR5 is also known as WD repeat-containing protein 5. These antibodies recognize WDR5 of human origin. Other species should be tested on a case-by-case basis.

### Control lysate information:

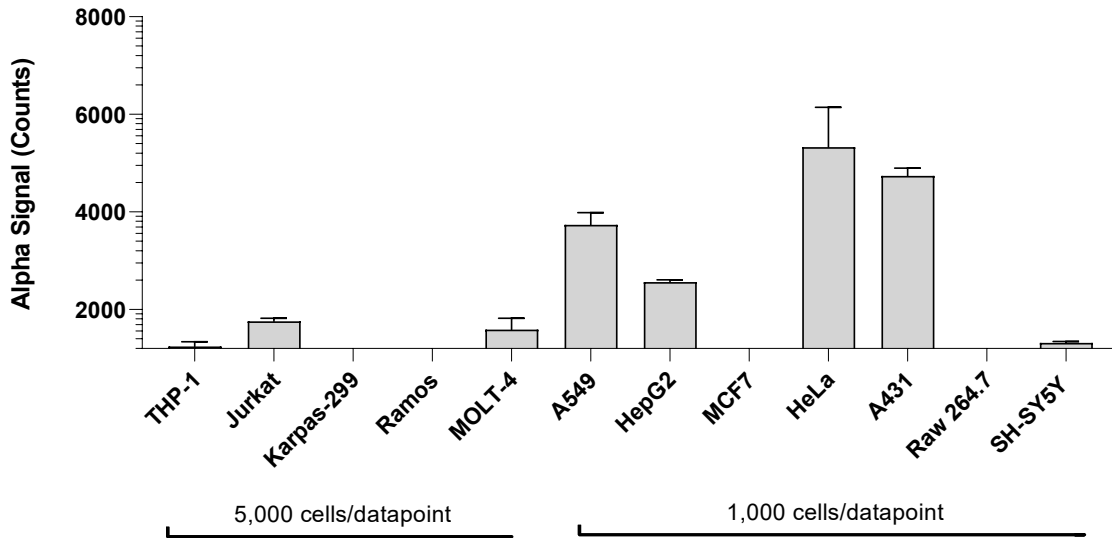
Positive Control Lysate: Prepared from Jurkat cells cultured in 10% FBS containing media. Cells were harvested and washed with HBSS + 0.1% BSA, adjusted to  $4 \times 10^6$ . Cells were then lysed with the addition of 5X Lysis Buffer B.

### Representative data

Data obtained with a 2-plate, 2-incubation protocol. A431 cells were plated at 20K cells/well in a 96 well plate and incubated overnight. Cells were treated with WDR5-IN-1 at the indicated concentrations for 20 hours. Cells were lysed with Lysis Buffer B (WDR5) or Lysis Buffer (Cofilin) and assayed separately for WDR5 Total and Cofilin Total using respective *SureFire Ultra* kits. Equivalent to approximately 2,000 cells/datapoint (WDR5 Total) or 40 cells/datapoint (Cofilin).



Data obtained from measurement of WDR5 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"

