

Research use only. Not for use in diagnostic procedures.

AlphaLISA[®] SureFire[®] Ultra[™]

Human and Mouse BRD9 Total Detection Kit

Product number:

ALSU-TBRD9-A500, ALSU-TBRD9-A10K,

ALSU-TBRD9-A50K, ALSU-TBRD9-A-HV



Kit specificity:

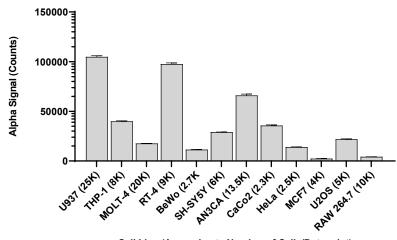
This assay kit contains antibodies which recognize distinct epitopes on BRD9. The protein detected by this kit corresponds to UniProt ID Q9H8M2. BRD9 is also known as Bromodomain-containing protein 9 and Rhabdomyosarcoma antigen MU-RMS-40.8. These antibodies recognize BRD9 of human and mouse origin. Other species should be tested on a case-by-case basis.

Control lysate information:

Positive Control Lysate: Prepared from U-937 cells, cultured to confluence in T175 flasks in 10% FBS containing media. Cells were harvested and washed in HBSS + 0.1% BSA, and lysed with Lysis Buffer at 7.5 x 10⁶ cells/mL.

Representative data

Data obtained from measurement of BRD9 Total in various cell types lysed with Lysis Buffer. Approximate number of cells/datapoint is indicated for the various cell lines.



Cell Line (Approximate Number of 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"





Revvity, Inc. 940 Winter Street Waltham, MA 02451 USA (800) 762-4000 www.revvity.com For a complete listing of our global offices, visit <u>www.revvity.com</u> Copyright ©2023, Revvity, Inc. All rights reserved.