

AlphaLISA® SureFire® Ultra™

Human and Mouse HDAC5 Total Detection Kit

Product number: ALSU-THDAC5-A500, ALSU-THDAC5-A10K,
ALSU-THDAC5-A50K, ALSU-THDAC5-A-HV



Kit specificity:

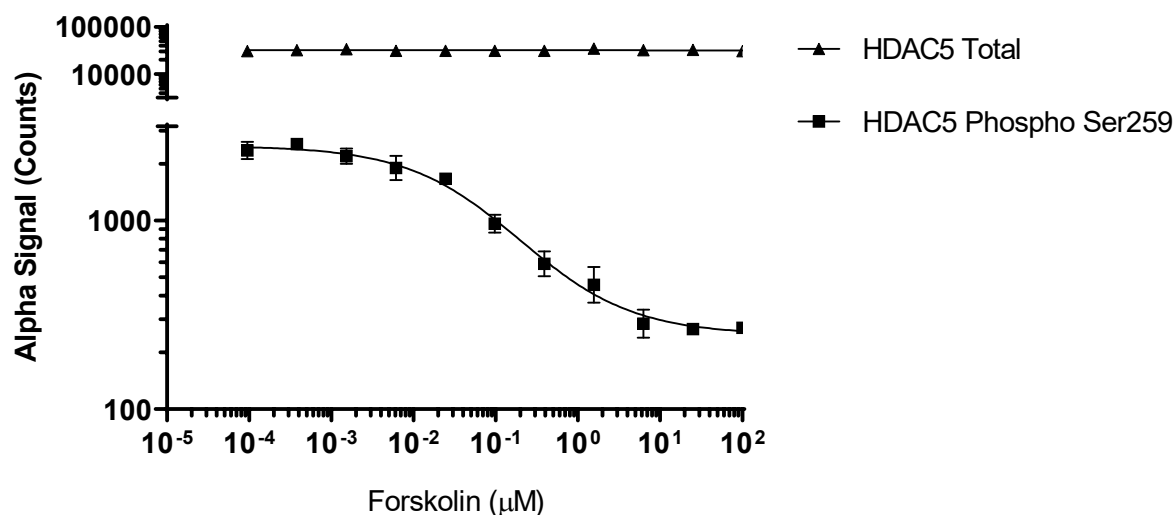
This assay kit contains antibodies which recognize distinct epitopes on HDAC5. The protein detected by this kit corresponds to UniProt ID Q9UQL6. HDAC5 is also known as histone deacetylase 5. These antibodies recognize HDAC5 of human and mouse origin. Other species should be tested on a case-by-case basis.

Control lysate information:

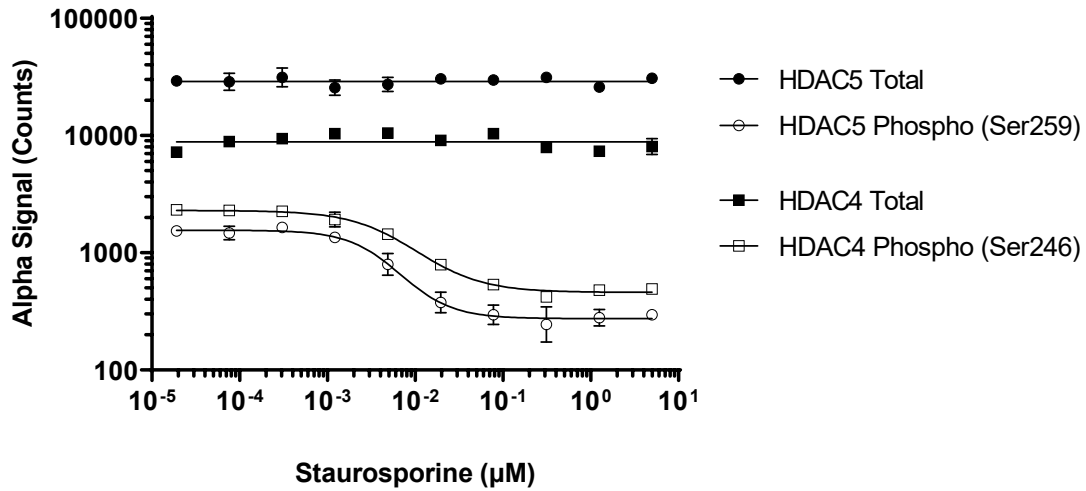
Positive Control Lysate: Prepared from NIH/3T3 cells, cultured to confluence in T175 flasks in 10% FBS containing media, then treated with 50 nM Calyculin A for 1 hour and lysed with 4 mL of Lysis Buffer.

Representative data

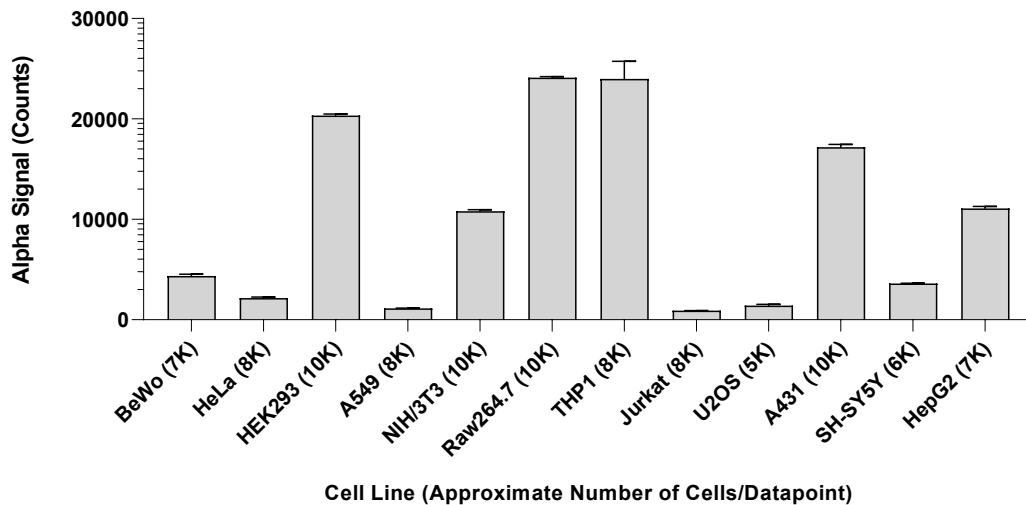
Data obtained with a 2-plate, 2-incubation protocol. HEK293 cells were plated at 20K cells/well in a 96 well plate and incubated for 48 hours. Cells were treated with Forskolin at the indicated concentrations for 1 hour. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Ser259) and Total HDAC5 using respective *SureFire Ultra* kit. Equivalent to approximately 8,000 cells/datapoint.



Data obtained with a 2-plate, 2-incubation protocol. NIH/3T3 cells were plated at 20K cells/well in a 96 well plate and incubated for 48 hours. Cells were treated with Staurosporine at the indicated concentrations for 1 hour. Cells were lysed with Lysis Buffer and assayed separately for various targets using respective *SureFire Ultra* kit. Equivalent to approximately 8,000 cells/datapoint.



Data obtained from measurement of HADC5 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"

