

# Technical Data Sheet

## AlphaLISA® SureFire® Ultra™ HDAC4 Total Assay Kit

ALSU-THDAC-A500, ALSU-THDAC-A10K, ALSU-THDAC-A50K, ALSU-THDAC-A-HV

### Kit Specificity

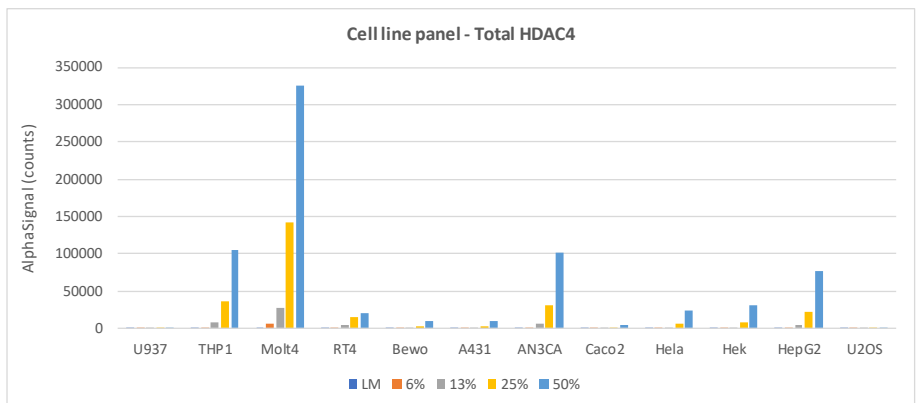
This assay kit contains antibodies which recognize invariant epitopes on HDAC4. The protein detected by this kit corresponds to UniProt ID P56524. HDAC4 is also known as Histone deacetylase 4, Histone deacetylase A, HD4 and HDACA. These antibodies recognize HDAC4 of human and mouse origin. Other species should be tested on a case-by-case basis.

### Control Lysate Information

Positive Control Lysate: Prepared from MOLT-4 cells cultured in T175 flasks to  $1.5 \times 10^6$  cells/mL in medium containing 10% FBS, and lysed in 1x *SureFire Ultra* Lysis Buffer at a final cell concentration of  $10 \times 10^6$  cells/mL.

### Representative Data

Data obtained with a 2-plate, 2-incubation protocol. A dilution series of various cell lysates from a panel of cell lines were prepared in 1X *SureFire Ultra* Lysis buffer at the approximate cell numbers indicated below. Lysates were assayed for Total HDAC4 by *SureFire Ultra*.



Cell Type	U937	THP1	Molt4	RT4	Bewo	A431	AN3CA	Caco2	HeLa	HEK	HEPG2	U2OS
Cells/datapoint at "50%"	6,000	16,000	40,000	18,000	5,500	15,000	27,500	4,500	16,000	20,000	14,000	10,000

For more information and detailed protocols, please refer to the generic AlphaLISA *SureFire Ultra* Assay Kits Manual, available from <http://www.perkinelmer.com/category/alpha-surefire-kits>.