

# Technical Data Sheet

## AlphaLISA® SureFire® Ultra™ Beclin-1 Total Assay Kit

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

### Kit Specificity

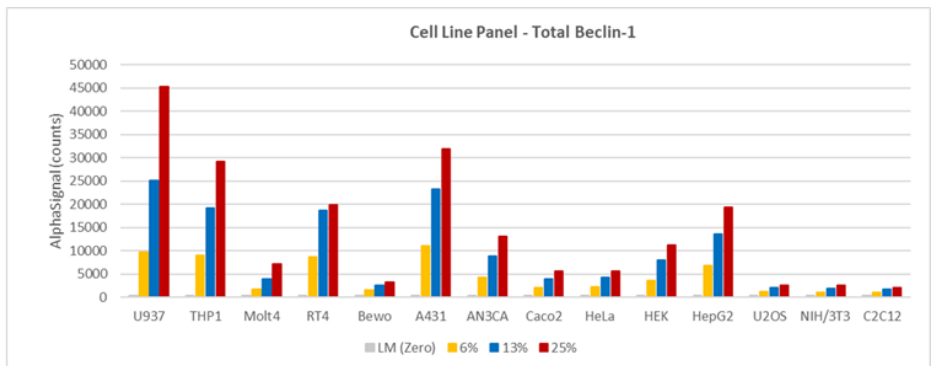
This assay kit contains antibodies which recognize invariant epitopes on Beclin-1. The protein detected by this kit corresponds to UniProt ID Q14457. Beclin-1 is also known as ATG-6. These antibodies recognize Beclin-1 of human, mouse and rat origin. Other species should be tested on a case-by-case basis.

### Control Lysate Information

Positive Control Lysate: Prepared from U937 cells, seeded at 800 K/mL and grown overnight in T175 flasks in 50 mL of 10% FBS containing medium. Cells were harvested, washed with HBSS and lysed in 5 mL of 1X *SureFire Ultra* Lysis Buffer.

### Representative Data

Data obtained for Total Beclin-1 from screening various lysates prepared from cells lysed in *SureFire Ultra* Lysis Buffer. The '25%' point corresponds to the approximate number of cells/datapoint indicated for the various cell lines.



Cell Type	U937	THP1	Molt4	RT4	Bewo	A431	AN3CA	Caco2	HeLa	HEK	HEPG2	U2OS	NIH3T3	C2C12
Cells/datapoint at "25%"	25,000	8,000	20,000	9,000	2,750	7,500	13,750	2,250	8,000	10,000	7,000	5,000	4,000	8,000

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>