

AlphaLISA® SureFire® Ultra™

## Human and Mouse p-ATG14 (Ser29) Detection Kit

**Product number:** ALSU-PATG14-A500, ALSU-PATG14-A10K,  
 ALSU-PATG14-A50K, ALSU-PATG14-A-HV



### Kit specificity:

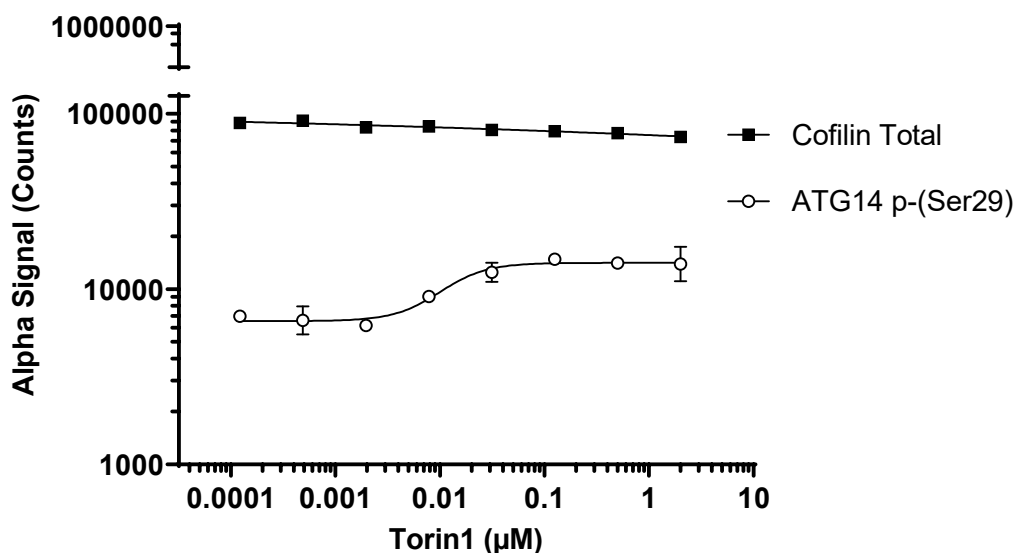
This assay kit contains antibodies which recognize the phospho Ser29 epitope and a distal epitope on ATG14. The protein detected by this kit corresponds to UniProt ID Q6ZNE5. ATG14 is also known as Beclin 1-associated autophagy-related key regulator and Barkor. These antibodies recognize ATG14 of human and mouse origin. Other species should be tested on a case-by-case basis.

### Control lysate information:

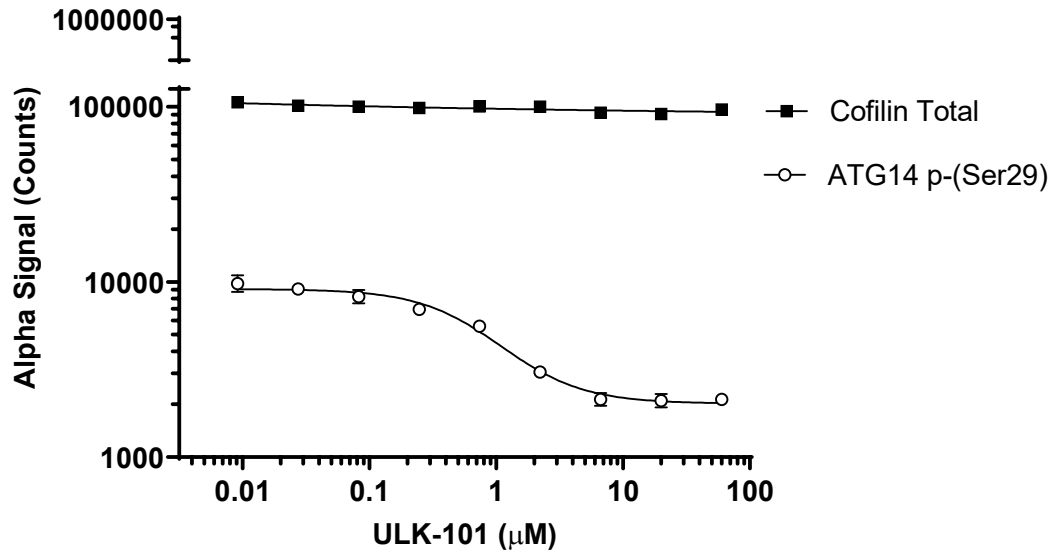
Positive Control Lysate: Prepared from PC-3 cells, cultured to confluence in T175 flasks in 10% FBS containing medium. Cells were treated with 200 nM Torin 1 for 3 hours and lysed with 2 mL Lysis Buffer.

### Representative data

Data obtained with a 2-plate, 2-incubation protocol. PC-3 cells were seeded at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with Torin 1 at the indicated concentrations for 3 hours. Cells were washed in HBSS, lysed with Lysis Buffer and assayed separately for Phospho ATG14 (Ser29) and Total Cofilin using respective *SureFire Ultra* kits. Equivalent to approximately 8,000 cells/datapoint (160 cells/datapoint for Cofilin Total).



Data obtained with a 2-plate, 2-incubation protocol. PC-3 cells were seeded at 40K cells/well in a 96 well plate and incubated overnight. Cells were then treated with ULK-101 at the indicated concentrations for 3 hours. Cells were washed with HBSS, lysed with Lysis Buffer and assayed separately for Phospho ATG14 (Ser29) and Total Cofilin using respective *SureFire Ultra* kits. Equivalent to approximately 8,000 cells/datapoint (160 cells/datapoint for Cofilin Total).



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"

