

AlphaLISA® SureFire® Ultra™

## Human and Mouse p-c-Myc (Ser62) Detection Kit

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



### Kit specificity:

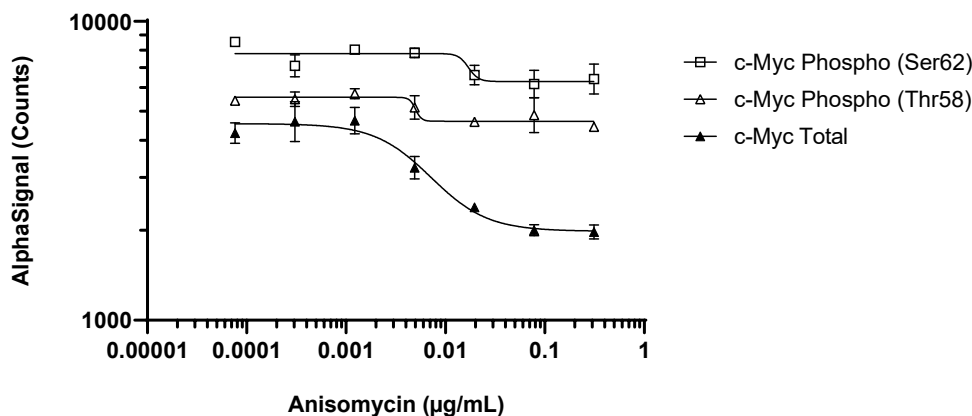
This assay kit contains antibodies which recognize the phospho-Ser62 epitope, and a distal epitope on c-Myc. The protein detected by this kit corresponds to UniProt ID P01106. c-Myc is also known as Transcription factor p64 and bHLHe39. These antibodies recognize c-Myc of human and mouse origin. Other species should be tested on a case-by-case basis.

### Control lysate information:

Positive Control Lysate: Prepared from HeLa cells, cultured to confluence in T175 flasks in 10% FBS containing media, treated with 10 $\mu$ M MG132 and 2 $\mu$ g/mL PMA for 30 minutes and lysed with 5 mL of Lysis Buffer.

### Representative data

Data obtained with a 2-plate, 2-incubation protocol. HeLa cells were seeded at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with Anisomycin at the indicated concentrations in the presence of 10  $\mu$ M MG-132 for 30 minutes. Cells were lysed with Lysis Buffer and assayed for various targets using respective *SureFire Ultra* kits. Equivalent to approximately 4000 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"