

# Technical Data Sheet

Caution: Research use only.

## AlphaLISA® SureFire® Ultra™ c-Myc Total Assay Kit

**Products: ALSU-TCMYC-A500, ALSU-TCMYC-A10K, ALSU-TCMYC-A50K,  
ALSU-TCMYC-A-HV**

### **Kit-Specificity Information**

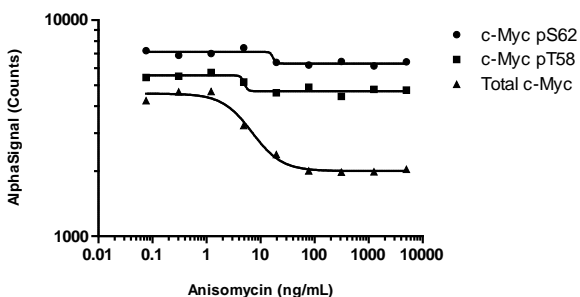
This assay kit contains antibodies which recognize invariant epitopes on c-Myc. The protein detected by this kit corresponds to Entrez Gene ID 4609. C-Myc is also known as avian myelocytomatosis viral oncogene homolog, BHLHE39 and bHLHe39MRTL. These antibodies recognize c-Myc of human origin. Other species 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 medium, then treated with 2 µg/mL PMA and 10 µM MG-132 for 30 min and lysed in 5 mL of 1X *SureFire Ultra* Lysis buffer.

### **Representative data**

Data obtained with a 2-plate, 2-incubation protocol. HeLa cells were plated overnight at 40K cells/well, then inhibited with the indicated concentrations of Anisomycin and 10 µM MG-132 for 30 min. Cells were lysed with 100 µL of 1X Lysis Buffer and assayed separately for total and phospho c-Myc, using respective kits.



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>