

# Technical Data Sheet

Caution: Research use only.

## AlphaLISA® SureFire® Ultra™ p-p53 (Ser15) Assay Kit

**Products: ALSU-PP53-A500, ALSU-PP53-A10K, ALSU-PP53-A50K,  
ALSU-PP53-A-HV**

### Kit-Specificity Information

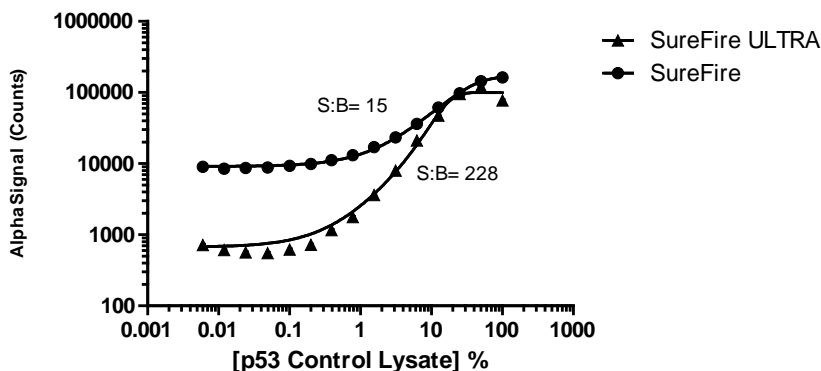
This assay kit contains antibodies which recognize the phospho-Ser15 epitope, and a distal epitope on p53. The protein detected by this kit corresponds to GenBank Accession NP\_000537. p53 is also known as P53, LFS1, TRP53, FLJ92943 and TP53. These antibodies recognize p53 of human origin. Other species should be tested on a case-by-case basis.

### Control Lysate Information

Positive Control Lysate: Prepared from U2OS cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with Doxorubicin for 18 hours, and lysed with 1X *SureFire Ultra* Lysis buffer.

### Representative data

Data obtained with a 2-plate, 2-incubation protocol. Control lysates, as indicated above, were serially diluted with 1X Lysis buffer from either *SureFire Ultra* or *SureFire* kits, and assayed with their respective kit. Data shows the much improved range and sensitivity of the *SureFire Ultra* assay of p53 (Ser15) phosphorylation compared to the *SureFire* assay.



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>