

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex p-p53 (Ser15) + Total p53 Assay Kit

Products: MPSU-PTP53-K500, MPSU-PTP53-K10K, MPSU-PTP53-K50K,
MPSU-PTP53-K-HV

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho Ser15 epitope, and a distal epitope, on p53, plus antibodies to measure the total levels of p53.

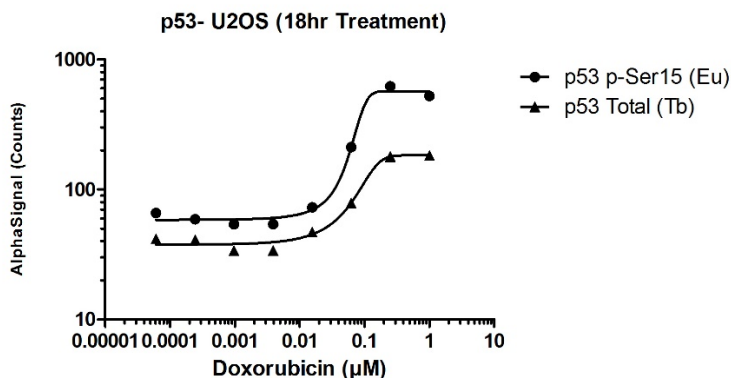
The protein detected by this kit corresponds to GenBank Accession NP_000537. p53 is also known as P53, LFS1, TRP53, FLJ92943 and TP53. These antibodies recognize p53 of human origin. Other species should be tested on a case-by-case basis.

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. Confluent U2OS cells were treated with Doxorubicin at the indicated concentrations for 18 hours, and lysed with 1X *SureFire Ultra* Lysis buffer. Samples were then assayed for both phospho p53 (Ser15) and total levels of p53 as detailed in the MPSU manual.



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>