

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ Ubiquitin Total Assay Kit

ALSU-TUBI-A500, ALSU TUBI-A10K, ALSU-TUBI-A50K, ALSU-TUBI-A-HV

Kit Specificity

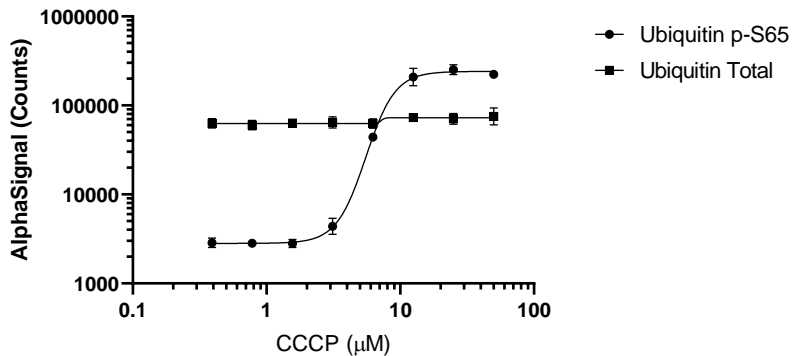
This assay kit contains antibodies which recognize invariant epitopes on Ubiquitin. The protein detected by this kit corresponds to UniProt ID P0CG47/P0CG48. Ubiquitin is also known as Ubiquitin B, Ubiquitin C, Polyubiquitin B, Polyubiquitin C, UBB & UBC. These antibodies recognize Ubiquitin of human and mouse origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from PC3 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 50 μ M carbonyl cyanide 3-chlorophenylhydrazone (CCCP) for 120 min and lysed in 4 mL of 1X *SureFire Ultra* Lysis Buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. PC3 cells were plated overnight at 30K cells/well, then stimulated with the indicated concentrations of carbonyl cyanide 3-chlorophenylhydrazone (CCCP) for 2 h. Cells were lysed with 100 μ L 1X *SureFire Ultra* Lysis Buffer and assayed separately for Total and Phospho (S65) Ubiquitin, using respective kits. Equivalent to approximately 3000 cells/datapoint.



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>