

AlphaLISA® SureFire® Ultra™

Human p-MDM2 (Ser166) Detection Kit

Product number: ALSU-PMDM2-A500, ALSU-PMDM2-A10K,
ALSU-PMDM2-A50K, ALSU-PMDM2-A-HV



Kit specificity:

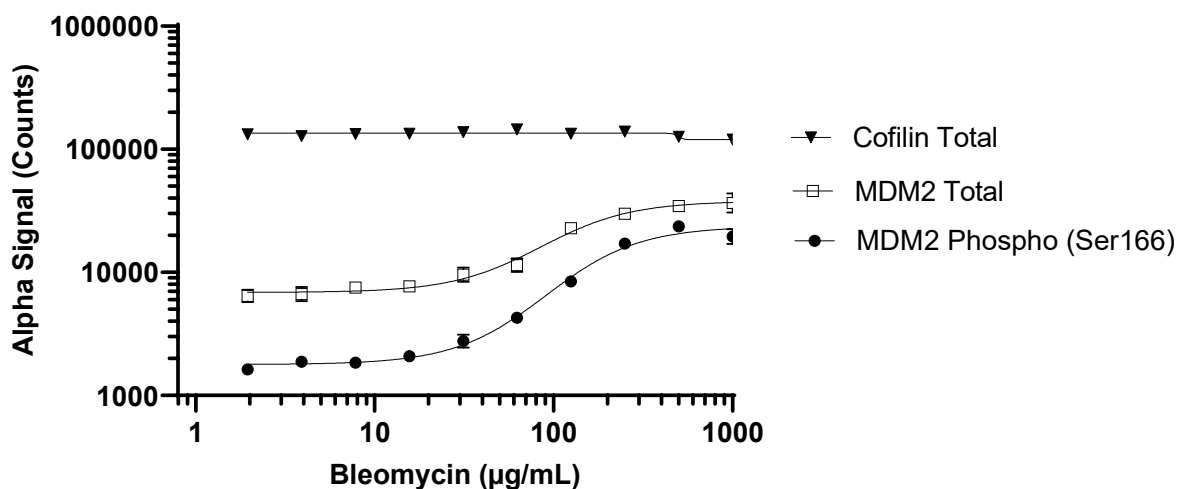
This assay kit contains antibodies which recognize the phospho-Ser166 epitope and a distal epitope on MDM2. The protein detected by this kit corresponds to UniProt ID F5H1M7. These antibodies recognize MDM2 of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

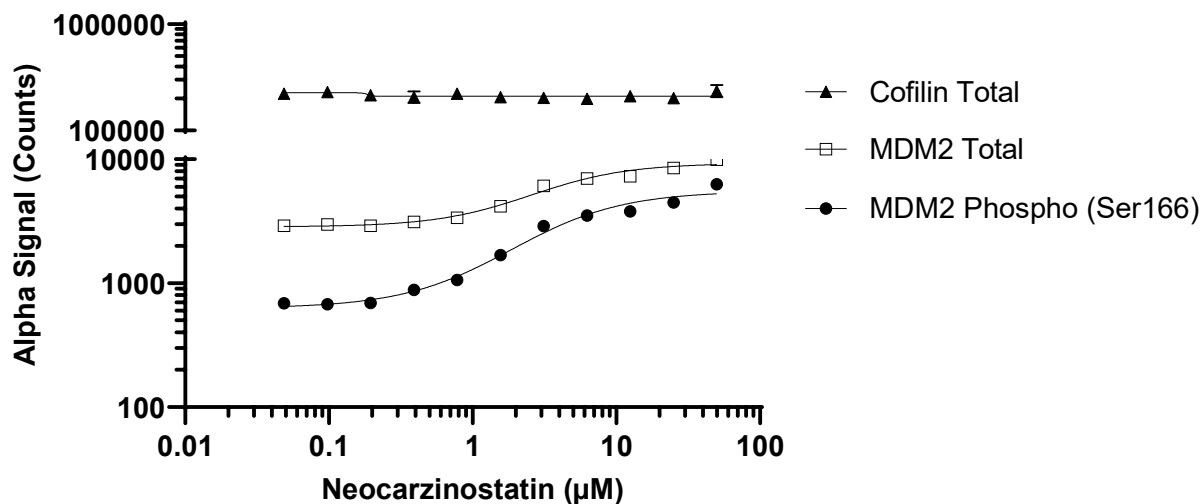
Positive Control Lysate: Prepared from MCF7 cells, cultured to confluence in T175 flasks in 10% FBS containing media, treated with 5 µM MG132 for 18 hours and lysed with 4 mL of Lysis Buffer B.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. MCF7 cells were seeded at 50K cells/well in a 96 well plate and incubated overnight. Cells were treated with Bleomycin at the indicated concentrations for 5 hours. Cells were lysed with Lysis Buffer B or Lysis Buffer for Cofilin Total and assayed separately for Phospho (Ser166) and Total MDM2, and Cofilin Total using respective *SureFire Ultra* kits. Equivalent to approximately 5,000 cells/datapoint or 500 cells/datapoint (Cofilin Total).



Data obtained with a 2-plate, 2-incubation protocol. MCF7 cells were seeded at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with Neocarzinostatin at the indicated concentrations for 7 hours. Cells were lysed with Lysis Buffer B or Lysis Buffer for Cofilin and assayed separately for Phospho (Ser166) and Total MDM2, and Cofilin Total using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint or 400 cells/datapoint (Cofilin Total).



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"

