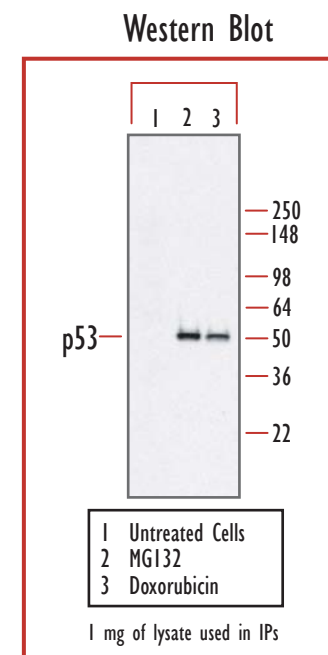
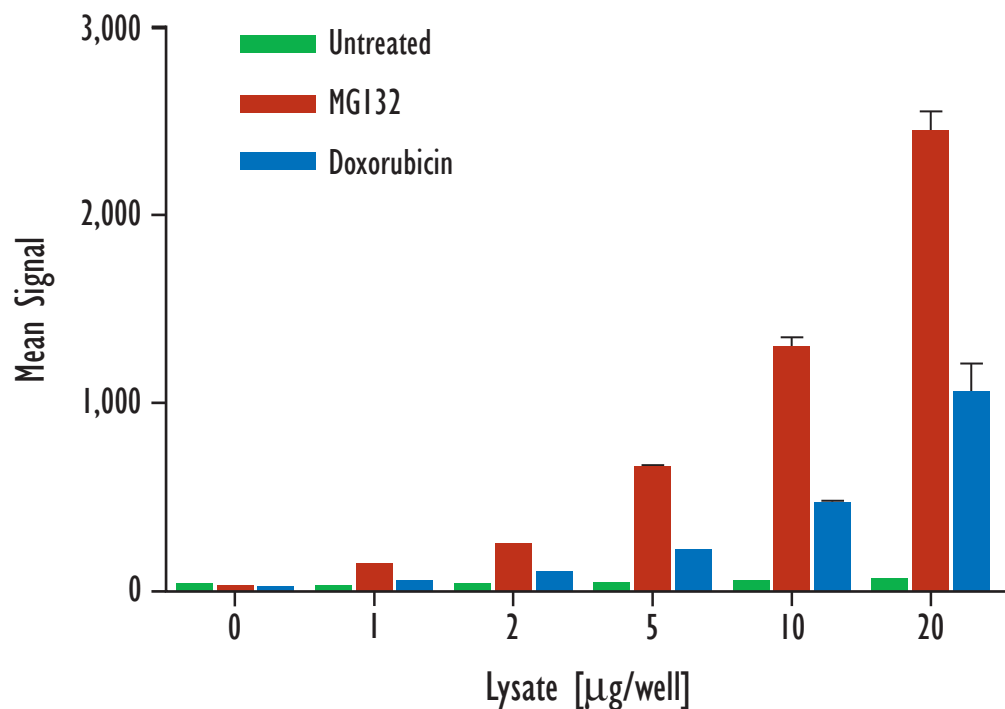
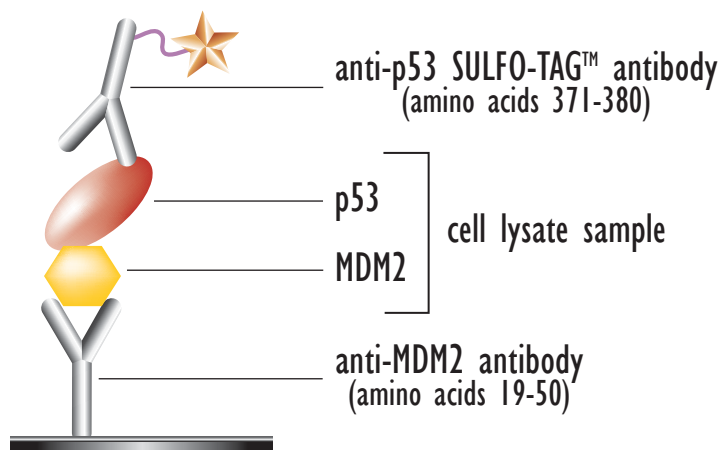


Detection of MDM2-p53 Complex



HCT116 cells were harvested after no treatment, or following treatment with MG132 or doxorubicin. Cells were lysed in MSD Complete Lysis Buffer. Whole cell lysates were diluted into MSD Complete Lysis Buffer and added to MSD MULTI-ARRAY™ plates coated with anti-MDM2 antibody. The MDM2-p53 complex was detected using anti-p53 antibody labeled with MSD SULFO-TAG reagent.

Note: RIPA buffer is not recommended as it may disrupt the MDM2-p53 complex.