

PathHunter P53 Translocation Assay

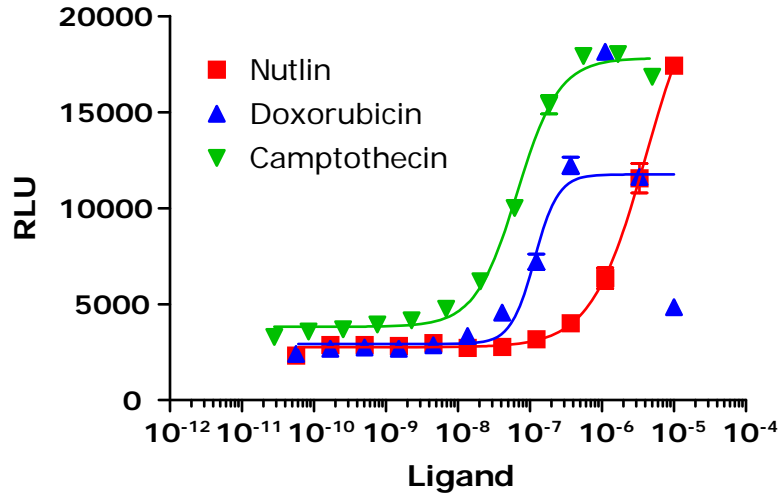


Figure 1. PathHunter® EA-Nuc parental cells were used to build P53 translocation assay. P53 was tagged to ProLabel and the data shows P53 translocation in response to compounds that affect DNA damage pathway.