

Technical Data Sheet

Alpha SureFire® Ultra™ Multiplex p-Rb (Ser780) + Total Rb Assay Kit

MPSU-PTRB-K500, MPSU-PTRB-K10K, MPSU-PTRB-K50K, MPSU-PTRB-K-HV

Kit-Specificity Information

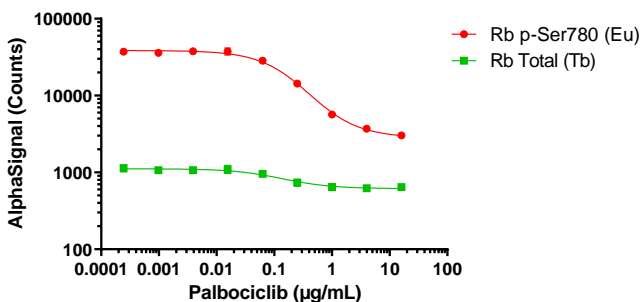
This assay kit contains antibodies which recognize the phospho-Ser780 epitope, and invariant epitopes on retinoblastoma tumor suppressor protein (Rb), plus antibodies to measure the total levels of Rb. The protein detected by this kit corresponds to UniProt ID P06400. Rb is also known as RB transcriptional corepressor 1. These antibodies recognize Rb of human origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from Jurkat cells cultured in T175 flasks to approximately 1×10^6 cells/mL in 10% FBS containing medium, then treated with 1.25 mM Pervanadate for 30 minutes and lysed at 3.2×10^6 cells/mL with 1X *SureFire Ultra* Lysis buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. THP-1 cells seeded into a 24 well plate at 2 million cells/mL and treated for 24 h with a dose range of Palbociclib. Cells were harvested, rinsed and lysed in *SureFire Ultra* Lysis Buffer. Lysates were screened for Phospho (Ser780) and Total Rb in the same wells of an assay plate by Alpha *SureFire Ultra* Multiplex. Equivalent to approximately 20,000 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>.