

AlphaLISA® SureFire® Ultra™ Multiplex

Human p-Rb (Ser807/811) + Total Rb Detection Kit

Product number: MPSU-PTRB-M500, MPSU-PTRB-M10K,
MPSU-PTRB-M50K, MPSU-PTRB-M-HV



Kit specificity:

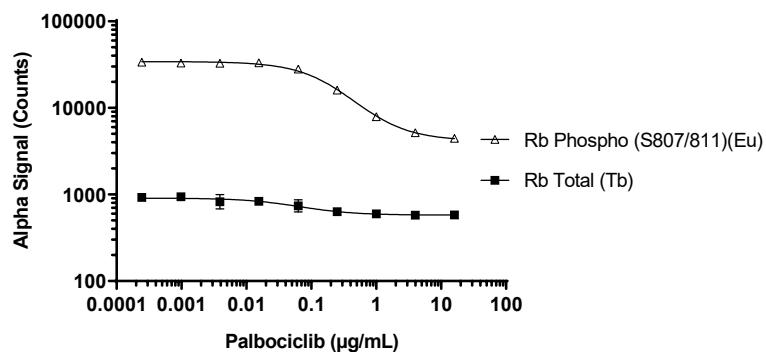
This assay kit contains antibodies which recognize the phospho-Ser807/811 epitope and a distal epitope on Rb, plus antibodies to detect Total Rb. The protein detected by this kit corresponds to UniProt ID P06400. Rb is also known as Retinoblastoma-associated protein and p110-RB1. These antibodies recognize 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 10% FBS containing medium. Cells were harvested and washed with HBSS + 0.1% BSA, adjusted to 4×10^6 cells/mL, then treated with 1.25 nM Pervanadate for 30 minutes. Cells were then lysed with 5X Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. THP-1 cells were seeded at 2×10^6 cells/mL in a 24 well plate and incubated for 20 minutes. Cells were treated with Palbociclib at the indicated concentrations for 24 hours. Cells were then washed with HBSS + 0.1% BSA, lysed with Lysis Buffer and assayed for Phospho (Ser807/811) and Total Rb in the same wells of an assay plate using the *SureFire Ultra* Multiplex kit. Equivalent to approximately 20,000 cells/datapoint.



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"