

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ Rb Total Assay Kit

**Products: ALSU-TRB-A500, ALSU-TRB-A10K, ALSU-TRB-A50K,
ALSU-TRB-A-HV**

Kit-Specificity Information

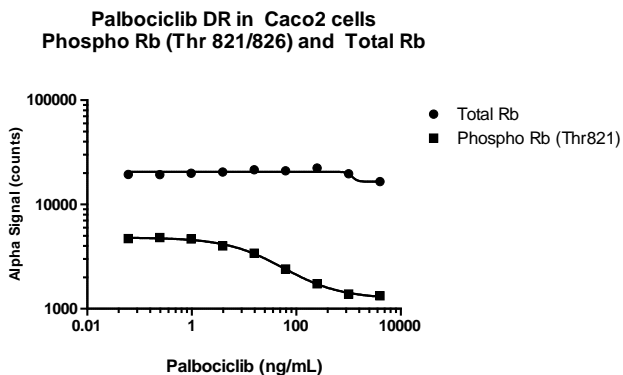
This assay kit contains antibodies which recognize invariant epitopes on retinoblastoma tumor suppressor protein (Rb). The human protein detected by this kit corresponds to Entrez Gene ID 5925. These antibodies recognize Rb of human origin. Other species should be tested on a case-by-case basis. Rb is also known as RB transcriptional corepressor 1.

Control Lysate Information

Positive Control Lysate: Prepared from Jurkat cells cultured in a T175 flask to 10^6 cells/mL in medium containing 10%FBS, treated with 2.5mM 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. Caco2 cells were seeded in a 96 well plate at 0.2×10^6 cells/mL, grown overnight, and treated a further 24 hours with Palbociclib at the indicated concentrations. Medium was removed and cells lysed with 100 μ L of 1X Lysis buffer. Samples were assayed separately for Total and Phospho (Thr821/826) Rb, using respective *SureFire Ultra* assay kits.



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>