Research use only. Not for use in diagnostic procedures.

AlphaLISA® SureFire® Ultra™

Human and Mouse p-p27 KIP1 (Ser10) Detection Kit

Product number: ALSU-PP27K-A500, ALSU-PP27K-A10K,

ALSU-PP27K-A50K, ALSU-PP27K-A-HV



Kit specificity:

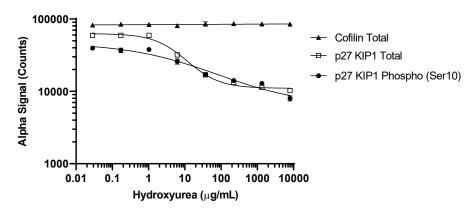
This assay kit contains antibodies which recognize the phospho-Ser10 epitope and a distal epitope on p27 KIP1. The protein detected by this kit corresponds to UniProt ID P46527. p27 KIP1 is also known as Cyclin-dependent kinase inhibitor 1B and CDKN1B. These antibodies recognize p27 KIP1 of human and mouse origin. Other species should be tested on a case-by-case basis.

Control lysate information:

Positive Control Lysate: Prepared from MCF7 cells, cultured to confluence in T175 flasks in 10% FBS containing media and lysed with 4 mL of Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. A431 cells were plated at 40K cells/well in a 96 well plate and incubated for 48 hours. Cells were treated with Hydroxyurea at the indicated concentrations for 18 hours. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Ser10), Total p27 KIP1 and Total Cofilin (housekeeping) using respective *SureFire Ultra* kits. Equivalent to approximately 8,000 cells/datapoint or 800 cells/datapoint for Total Cofilin.



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"



