

AlphaLISA® SureFire® Ultra™

Human p16 INK4A Total Detection Kit

Product number: ALSU-TP16INK4-A500, ALSU-TP16INK4-A10K,
 ALSU-TP16INK4-A50K, ALSU-TP16INK4-A-HV



Kit specificity:

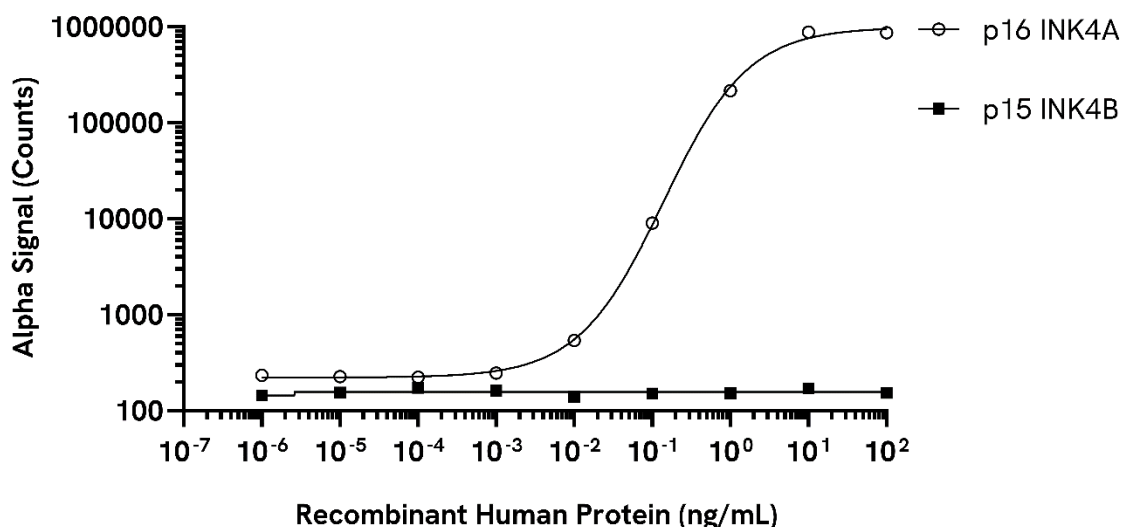
This assay kit contains antibodies which recognize distinct epitopes on p16 INK4A. The protein detected by this kit corresponds to UniProt ID P42771. p16 INK4A is also known as p16 INK4a, Cyclin-dependent kinase inhibitor 2A (CDKN2A) and Multiple tumor suppressor 1. These antibodies recognize p16 INK4A of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

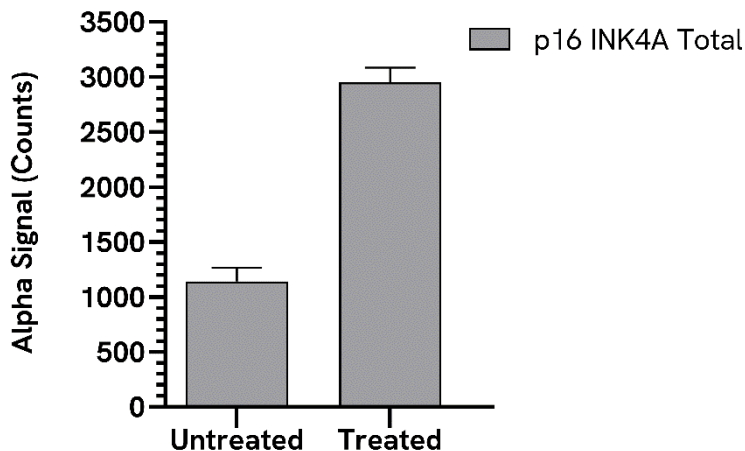
Positive Control Lysate: Prepared from HEK293 cells cultured to confluence in T175 flasks in 10% FBS containing medium and lysed with 8 mL of Lysis Buffer.

Representative data:

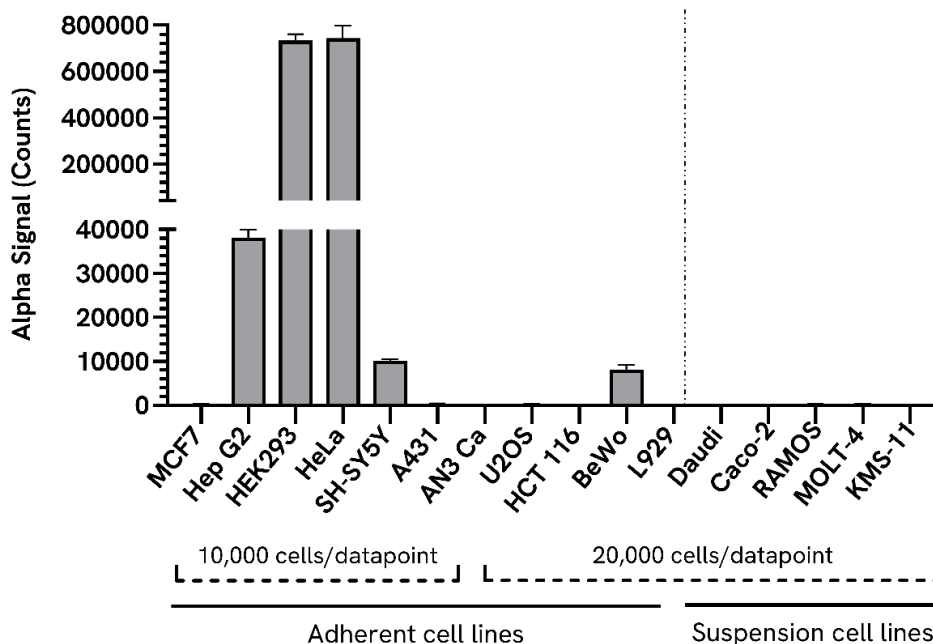
p16 INK4A (Abcam, Ab84075) and p15 INK4B (Abcam, ab268830) recombinant human proteins were serially diluted with Lysis Buffer and evaluated using the p16 INK4A Total SureFire Ultra kit. No cross-reactivity against p15 INK4B was observed despite sharing over 80% similarity with p16 INK4A.



Data obtained with a 2-plate, 2-incubation protocol. HEK293 cells were seeded at 10K cells/well in a 96 well plate and incubated overnight. Cells were treated with 2.5 μ M Palbociclib for 24 hours. Cells were lysed with Lysis Buffer and assayed for Total p16 INK4A using the *SureFire Ultra* kit. Equivalent to approximately 200 cells/datapoint.



Data obtained from measurement of p16 INK4A Total in various cell types lysed with Lysis Buffer. Approximate number of cells/datapoint is indicated for the various cell lines.



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"

