

AlphaLISA® SureFire® Ultra™

## Human and Mouse CDK5 Total Detection Kit

**Product number:** ALSU-TCDK5-A500, ALSU-TCDK5-A10K,  
ALSU-TCDK5-A50K, ALSU-TCDK5-A-HV



### Kit specificity:

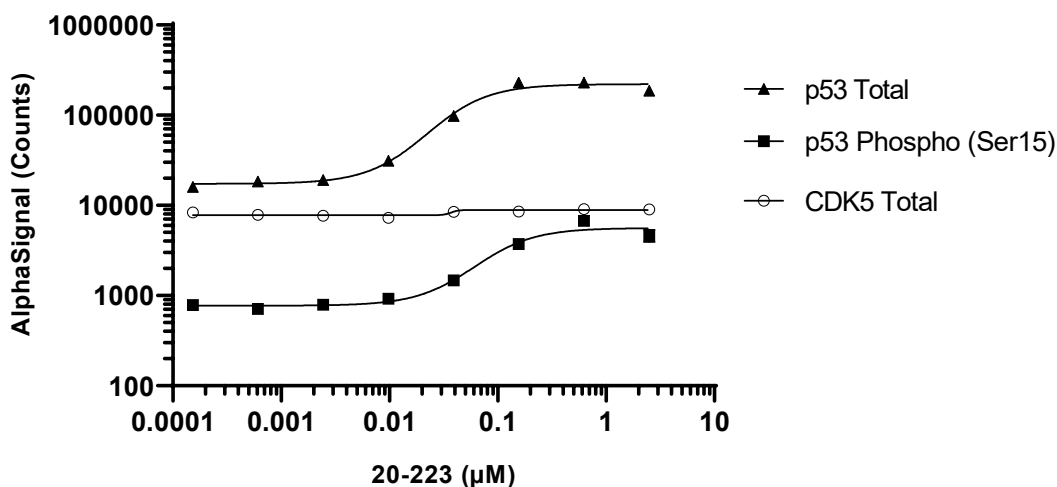
This assay kit contains antibodies which recognize distinct epitopes on CDK5. The protein detected by this kit corresponds to UniProt ID Q00535. CDK5 is also known as cyclin-dependent-like kinase 5. These antibodies recognize CDK5 of human and mouse origin. Other species should be tested on a case-by-case basis.

### Control lysate information:

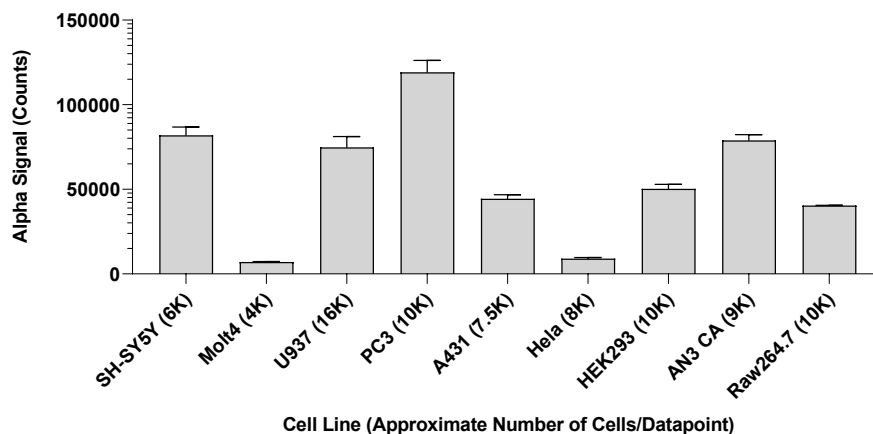
Positive Control Lysate: Prepared from HeLa cells, cultured to confluence in T175 flasks in 10% FBS containing media and lysed with 5 mL of Lysis Buffer.

### Representative data

Data obtained with a 2-plate, 2-incubation protocol. MCF7 cells were plated at 40K cells per well in a 96 well plate and incubated overnight. Cells were treated with CDK5 Inhibitor 20-223 at the indicated concentrations for 6 hours in the presence of 5 µg/mL Insulin. Cells were lysed with Lysis Buffer and assayed separately for Total CDK5 and Total and Phospho (Ser15) p53 using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint.



Data obtained from measurement of CDK5 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"

