

AlphaLISA® SureFire® Ultra™

## Human CDK6 Total Detection Kit

**Product number:** ALSU-TCDK6-A500, ALSU-TCDK6-A10K,  
ALSU-TCDK6-A50K, ALSU-TCDK6-A-HV



### Kit specificity:

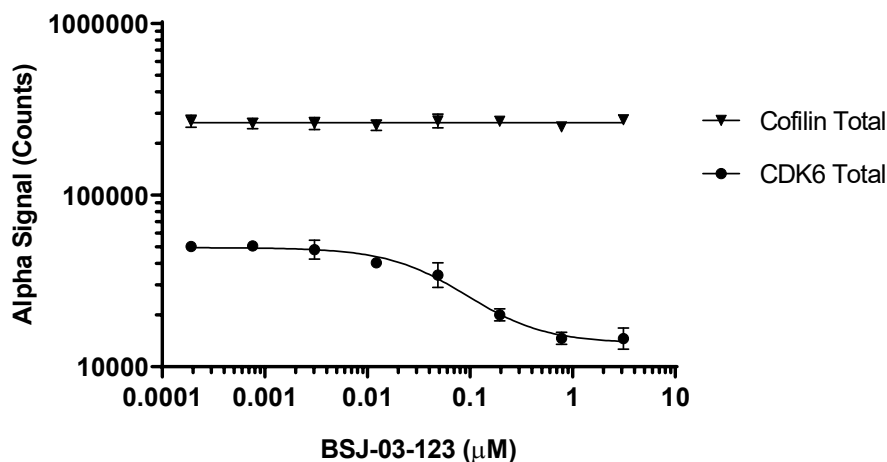
This assay kit contains antibodies which recognize distinct epitopes on CDK6. The protein detected by this kit corresponds to UniProt ID Q00534. CDK6 is also known as cyclin-dependent-like kinase 6. These antibodies recognize CDK6 of human origin. Other species should be tested on a case-by-case basis.

### Control lysate information:

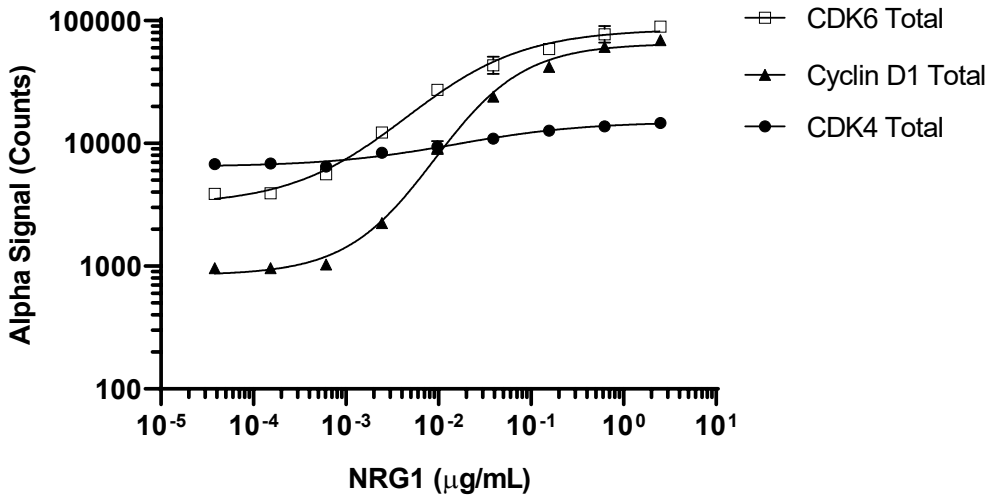
Positive Control Lysate: Prepared from MOLT-4 cells cultured in 10% FBS containing media. Cells were washed in HBSS + 0.1% BSA, spun down at 300 RCF for 5 minutes and then lysed in Lysis Buffer at  $2 \times 10^6$  cells/mL.

### Representative data

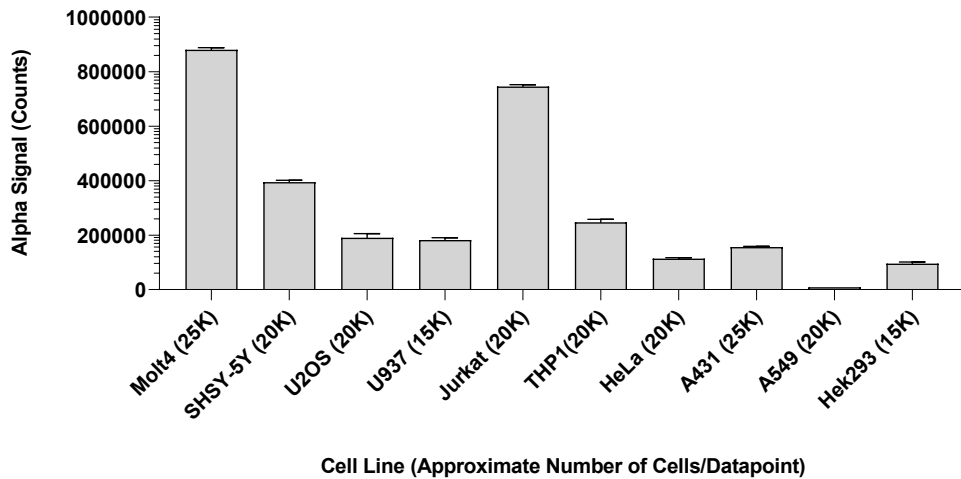
Data obtained with a 2-plate, 2-incubation protocol. SH-SY5Y cells were plated at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with BSJ-03-123 (CDK6 PROTAC Degradator) at the indicated concentrations for 6 hours. Cells were lysed with Lysis Buffer and assayed separately for Total CDK6 and Total Cofilin using respective *SureFire Ultra* kits. Equivalent to approximately 1,600 cells/datapoint.



Data obtained with a 2-plate, 2-incubation protocol. MCF7 cells were seeded at 20K cells/well in a 96 well plate and incubated overnight. Cells were treated with 20  $\mu$ M Wortmannin for 3 hours and further stimulated with Neuregulin1 at the indicated concentrations for 24 hours. Both treatments were performed in serum free media. Cells were lysed with Lysis Buffer and assayed separately for various targets using respective *SureFire Ultra* kits. Equivalent to approximately 2,000 cells/datapoint.



Data obtained from measurement of CDK6 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"

