

Research use only. Not for use in diagnostic procedures.

AlphaLISA™ SureFire® Biotin Free

Human p-CDK1 (Thr14) Detection Kit

Product number: ASBF-PCDK1-A500, ASBF-PCDK1-A10K,

ASBF-PCDK1-A50K, ASBF-PCDK1-A-HV



Kit specificity:

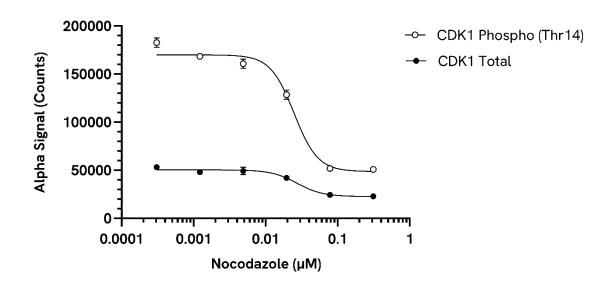
This assay kit contains antibodies which recognize the phospho-Thr14 epitope and a distal epitope on Cyclin-dependent kinase 1 (CDK1). The protein detected by this kit corresponds to UniProt ID P06493. CDK1 is also known as Cell division control protein 2 homolog and p34 protein kinase. These antibodies recognize CDK1 of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

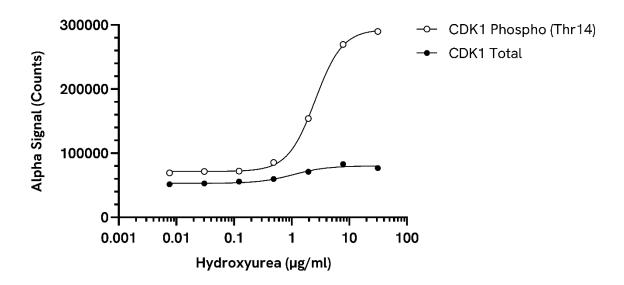
Positive Control Lysate: Prepared from HEK293 cells, cultured to confluence in T175 flasks in 10% FBS containing medium and lysed with 10 mL of Lysis Buffer.

Representative data:

Data obtained with a 2-plate, 2-incubation protocol. Jurkat cells were seeded at 200K cells/well in a 96-well plate in complete RPMI 1640 medium and treated with Nocodazole at the indicated concentrations for 18 hours. Cells were lysed with the addition of 5X Lysis Buffer and assayed separately for Phospho (Thr14) and Total CDK1 using respective *SureFire* Biotin Free kits. Equivalent to approximately 2,000 cells/datapoint.



Data obtained with a 2-plate, 2-incubation protocol. MOLT-4 cells were seeded at 200K cells/well in a 96-well plate in complete RPMI 1640 medium and treated with Hydroxyurea at the indicated concentrations for 18 hours. Cells were lysed with the addition of 5X Lysis Buffer and assayed separately for Phospho (Thr14) and Total CDK1 using respective *SureFire* Biotin Free kits. Equivalent to approximately 2,000 cells/datapoint.



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"



