

AlphaLISA® SureFire® Ultra™

Human DNA PKcs Total Detection Kit

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



Kit specificity:

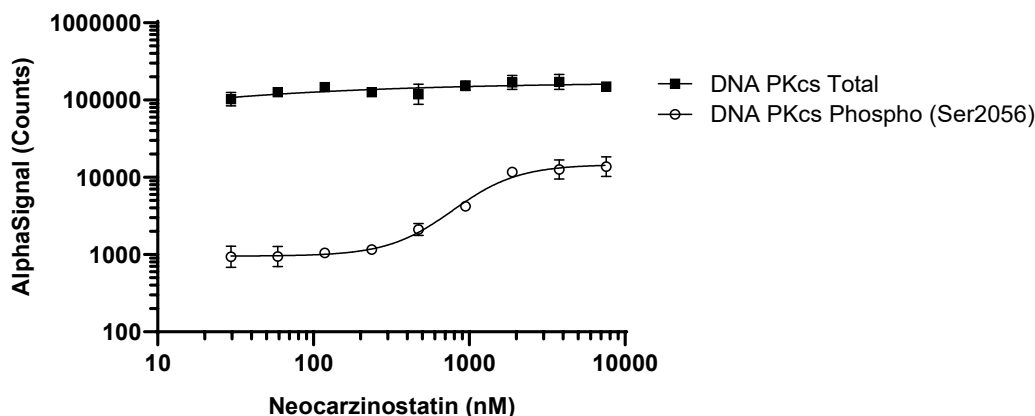
This assay kit contains antibodies which recognize distinct epitopes on DNA PKcs. The protein detected by this kit corresponds to UniProt ID P78527. DNA PKcs is also known as DNA-dependent protein kinase catalytic subunit, DNPk1 or p460. These antibodies recognize DNA PKcs of human origin. Other species 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, treated with 3 μM Neocarzinostatin for 2 hours and lysed with 4 mL of Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. RT4 cells were seeded at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with Neocarzinostatin at the indicated concentrations for 2 hours. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Ser2056) and Total DNA PKcs, using respective SureFire Ultra kits. Equivalent to approximately 4,000 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"