

Research use only. Not for use in diagnostic procedures.

AlphaLISA[®] SureFire[®] Ultra[™]

Human and Mouse KAP1 Total Detection Kit

Product number:

ALSU-TKAP1-A500, ALSU-TKAP1-A10K,

ALSU-TKAP1-A50K, ALSU-TKAP1-A-HV



Kit specificity:

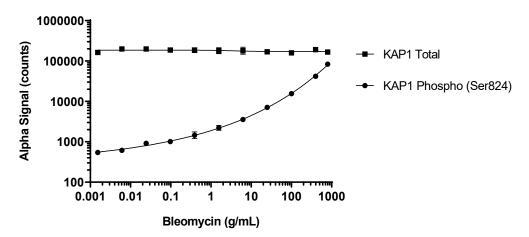
This assay kit contains antibodies which recognize distinct epitopes on KAP1. The protein detected by this kit corresponds to UniProt ID Q13263. KAP1 is also known as TIF1-beta, KRIP-1 and TRIM28. These antibodies recognize KAP1 of human and mouse origin. Other species should be tested on a case-by-case basis.

Control lysate information:

Positive Control Lysate: Prepared from HEK293 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, treated with 50 μ g/mL Bleomycin for 4 hours, and lysed in 8 mL Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. HeLa cells were plated at 20K cells per well and incubated overnight. Cells were then treated with Bleomycin at the indicated concentrations for 2 hours. Cells were lysed with Lysis Buffer and assayed separately for Phospho (S824) and Total KAP1 using respective *SureFire Ultra* 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"



Revvity, Inc. 940 Winter Street Waltham, MA 02451 USA (800) 762-4000 www.revvity.com For a complete listing of our global offices, visit <u>www.revvity.com</u> Copyright ©2023, Revvity, Inc. All rights reserved.

VersionTGRSU322.2

BioSciences