

Research use only. Not for use in diagnostic procedures.

AlphaLISA® SureFire® Ultra™

Human YAP Total Detection Kit

Product number: ALSU-TYAP-A500, ALSU-TYAP-A10K,

ALSU-TYAP-A50K, ALSU-TYAP-A-HV



Kit specificity:

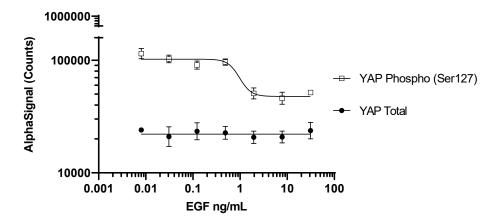
This assay kit contains antibodies which recognize distinct epitopes on YAP1. The protein detected by this kit corresponds to UniProt ID P46937. YAP is also known as Transcriptional coactivator YAP1 and Yes-associated protein 1. These antibodies recognize YAP1 of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

Positive Control Lysate: Prepared from AN3 CA 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. HeLa cells were plated at 40K cells/well in a 96 well plate and incubated overnight. Cells were serum starved overnight and then treated with EGF at the indicated concentrations for 60 minutes. Cells were lysed with Lysis Buffer and assayed for Phospho (Ser127) and Total YAP using respective SureFire Ultra assay 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"



