

# Technical Data Sheet

## AlphaLISA® SureFire® Ultra™ p-YAP1 (Ser127) Assay Kit

ALSU-PYAP-B500, ALSU-PYAP-B10K, ALSU-PYAP-B50K, ALSU-PYAP-B-HV

### Kit Specificity

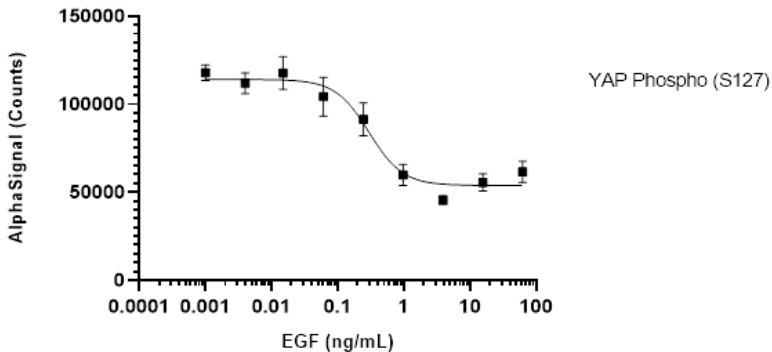
This assay kit contains antibodies which recognize the phospho Ser-127 epitope and a distal epitope on YAP1. The protein detected by this kit corresponds to UniProt ID P46937. YAP is also known as Transcriptional coactivator YAP1 or Yes-associated protein 1. These antibodies recognize YAP1 of mouse, rat and human origin. Other species should be tested on a case-by-case basis.

### Control Lysate Information

Positive Control Lysate: Prepared from AN3CA cells, cultured to confluence in T175 flasks in 10% FBS containing medium and lysed in 10 mL of 1X *SureFire Ultra* Lysis Buffer.

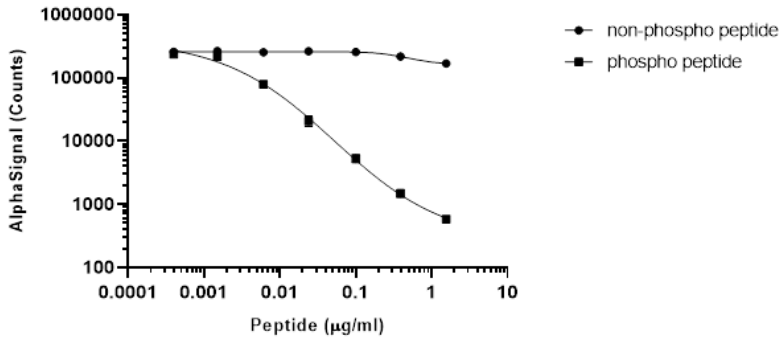
### Representative Data

Data obtained with a 2-plate, 2-incubation protocol. HeLa cells were plated at 20K cells per well in a 96 well plate and incubated overnight. Cells were serum starved overnight prior to stimulation with EGF at the indicated concentrations for 60 minutes. Cells were lysed with 1X *SureFire Ultra* Lysis Buffer and assayed for Phospho (Ser127) YAP using the *SureFire Ultra* assay kit. Equivalent to approximately 2000 cells/datapoint.



# Technical Data Sheet

Specificity of the assay was determined using peptides corresponding to the Ser127 epitope of YAP. Phosphorylated and non-phosphorylated peptides were titrated into control lysate and assayed for Phospho (Ser127) YAP, using the *SureFire Ultra* assay kit.



For more information and detailed protocols, please refer to the generic AlphaLISA *SureFire Ultra* Assay Kits Manual, available from <http://www.perkinelmer.com/category/alpha-surefire-kits>