

AlphaLISA® SureFire® Ultra™

Human AMPK α 1/2 Total Detection Kit

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



Kit specificity:

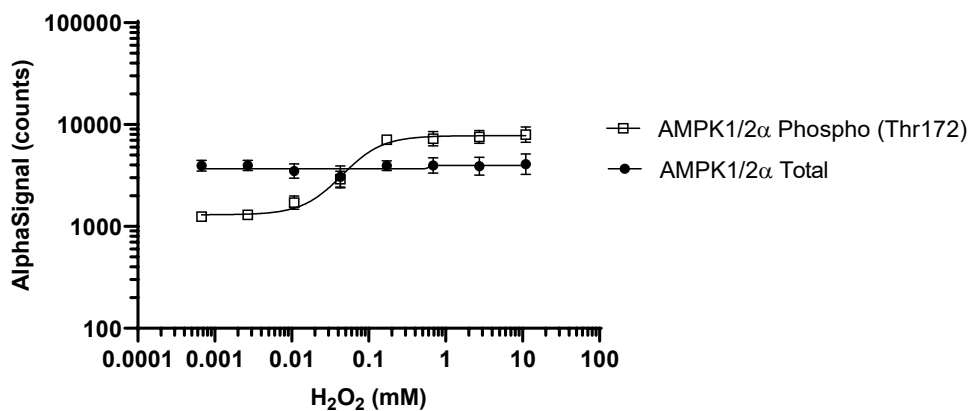
This assay kit contains antibodies which recognize distinct epitopes on AMPK α 1/2. The protein detected by this kit corresponds to UniProt ID Q13131 (AMPK α 1) and P54646 (AMPK α 2). AMPK α 1/2 is also known as AMP-activated protein kinase. These antibodies recognize AMPK α 1/2 of human origin. Other 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 media, treated with 11 mM H₂O₂ for 10 minutes and lysed with 5 mL of Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. HEK293 were seeded at 40K cells per well in a 96 well plate and incubated overnight. Cells were treated with H₂O₂ at the indicated concentrations for 10 minutes. Cells were lysed with Lysis Buffer and assayed for Phospho (Thr172) and Total AMPK α 1/2 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"