

Research use only. Not for use in diagnostic procedures.

AlphaLISA® SureFire® Ultra™ Human ACC Total Detection Kit

Product number:

ALSU-TACC-A500, ALSU-TACC-A10K,

ALSU-TACC-A50K, ALSU-TACC-A-HV



Kit specificity:

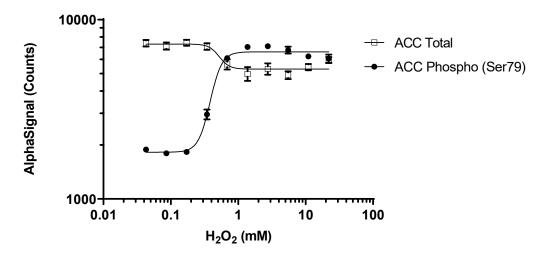
This assay kit contains antibodies that recognize distinct epitopes on ACC. The protein detected by this kit corresponds to UniProt ID Q13085. ACC is also known as Acetyl-coA carboxylase 1 (alpha). These antibodies recognize ACC of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

Positive control lysate: Prepared from HEK293 cells, cultured to confluence in T175 flasks in 10% FBS containing media, then 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 cells were seeded at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with H_2O_2 at the indicated concentrations for 10 minutes. Cells were lysed with Lysis Buffer and diluted further prior to assaying for Phospho (Ser79) and Total ACC, using respective *SureFire Ultra* assay kits. Equivalent to approximately 200 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.

VersionTGRSU290.2