

AlphaLISA® SureFire® Ultra™

Human CD3 Zeta Total Detection Kit

Product number: ALSU-TCD3-A500, ALSU-TCD3-A10K,
ALSU-TCD3-A50K, ALSU-TCD3-A-HV



Kit specificity:

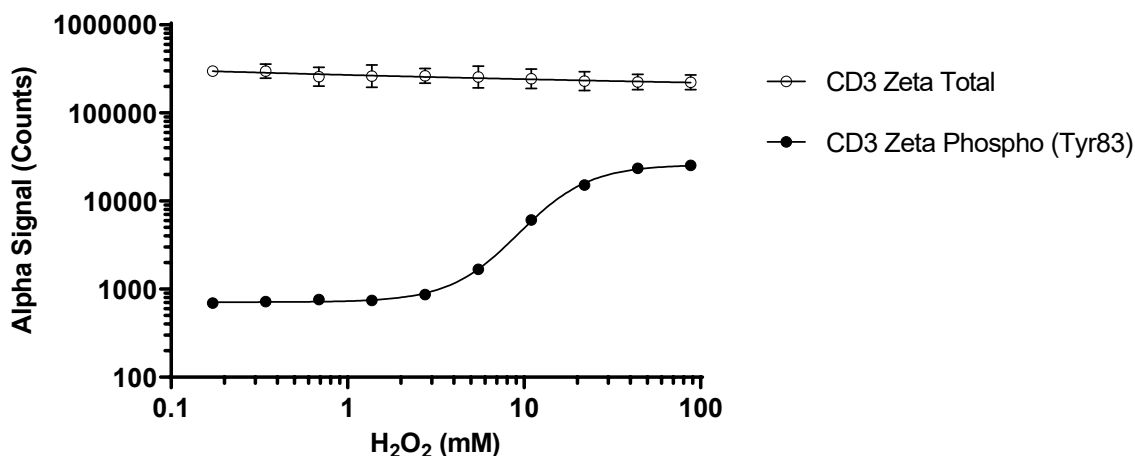
This assay kit contains antibodies which recognize distinct epitopes on CD3 Zeta. The protein detected by this kit corresponds to UniProt ID P20963. CD3 Zeta is also known as CD247. These antibodies recognize CD3 Zeta of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

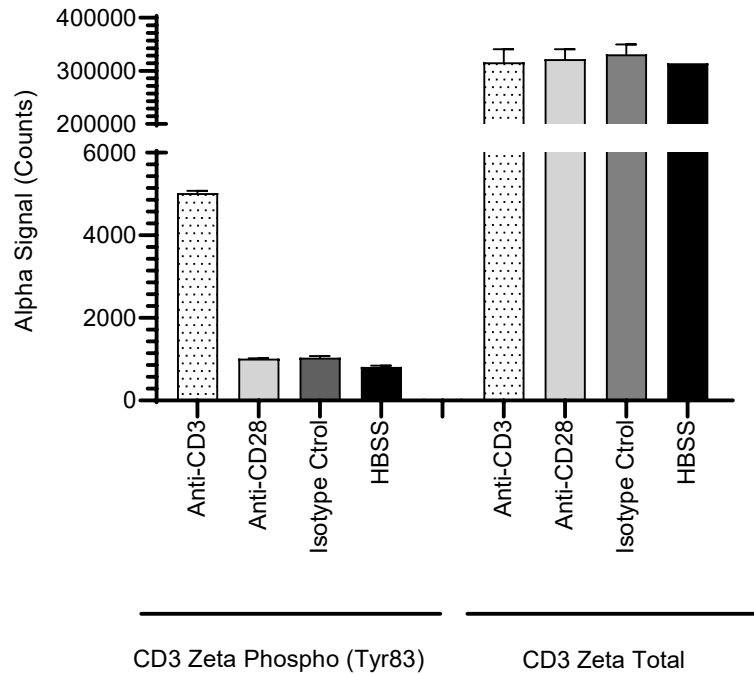
Positive Control Lysate: Prepared from Jurkat cells cultured in 10% FBS containing media. Cells were harvested and washed with HBSS + 0.1% BSA. Cells were adjusted to 3.5x10⁶ cells/mL and treated with 44 mM H₂O₂ for 10 minutes. Cells were then lysed with the addition of 5X Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. Jurkat cells were plated at 200K cells/well in a 96 well plate in HBSS + 0.1% BSA, then treated with H₂O₂ at the indicated concentrations for 10 minutes at 37°C. Cells were lysed with 5X Lysis Buffer and assayed separately for Phospho (Tyr83) and Total CD3 Zeta using respective SureFire Ultra kits. Equivalent to approximately 8,000 cells/datapoint.



Data obtained with a 2-plate, 2-incubation protocol. Jurkat cells were plated at 5×10^6 cells/mL in 12 well plates in HBSS + 0.1% BSA. Cells were treated for 5 minutes with $10 \mu\text{g/mL}$ of anti-CD3 Ab (BD cat# 555329) or a mouse IgG isotype control. Cells were lysed with the addition of 5X Lysis Buffer and lysates were further diluted in Lysis Buffer before assayed for Phospho (Tyr83) and Total CD3 Zeta. Equivalent to approximately 8,000 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"

