

AlphaLISA® SureFire® Ultra™

Human p-Lck (Tyr505) Detection Kit

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



Kit specificity:

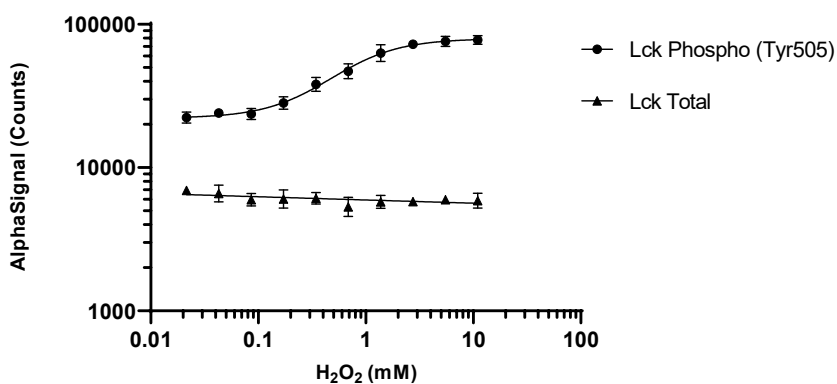
This assay kit contains antibodies which recognize the phospho-Tyr505 epitope and a distal epitope on Lck. The protein detected by this kit corresponds to UniProt ID P06239. Lck is also known as LSK and Protein YT16. These antibodies recognize Lck of human origin. Other species should be tested on a 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, adjusted to 4×10^6 cells/mL and then treated with 11 mM H₂O₂ and 50 nM Calyculin A 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 harvested, washed in HBSS and plated at a density of 1×10^6 cells/mL in a 96 well plate and then treated with H₂O₂ at the indicated concentrations for 10 minutes. Cells were lysed with 5X Lysis Buffer at a final cell density of 400K cells/mL. Lysates were further diluted and assayed for Phospho (Y505) and Total Lck using respective *SureFire Ultra* kits. Equivalent to approximately 800 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"