

AlphaLISA® SureFire® Ultra™

## Human p-PKCθ (Thr538) Detection Kit

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



### Kit specificity:

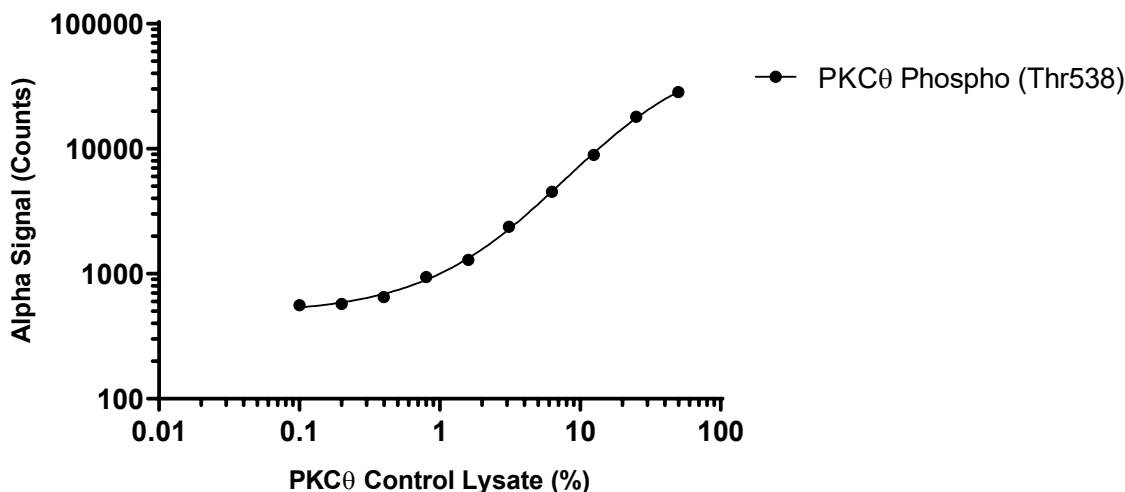
This assay kit contains antibodies which recognize the phospho-Thr538 epitope and a distal epitope on PKC-theta (PKCθ). The protein detected by this kit corresponds to UniProt ID Q04759. These antibodies recognize PKCθ of human origin. Other species should be tested on a case-by-case basis.

### Control lysate information:

Positive Control Lysate: Prepared from Jurkat cells cultured in 10% FBS containing media. Cells were washed with HBSS + 0.1% BSA, then spun down at 300 RCF for 5 minutes and lysed in Lysis Buffer at  $4 \times 10^6$  cells/mL.

### Representative data

Data obtained with a 2-plate, 2 incubation protocol. Jurkat control lysate was diluted to the indicated concentrations and assayed for Phospho (Thr538) PKCθ using the *SureFire Ultra* kit.



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"