

AlphaLISA® SureFire® Ultra™

Human and Mouse p-CaMKII (Thr286) Detection Kit

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



Kit specificity:

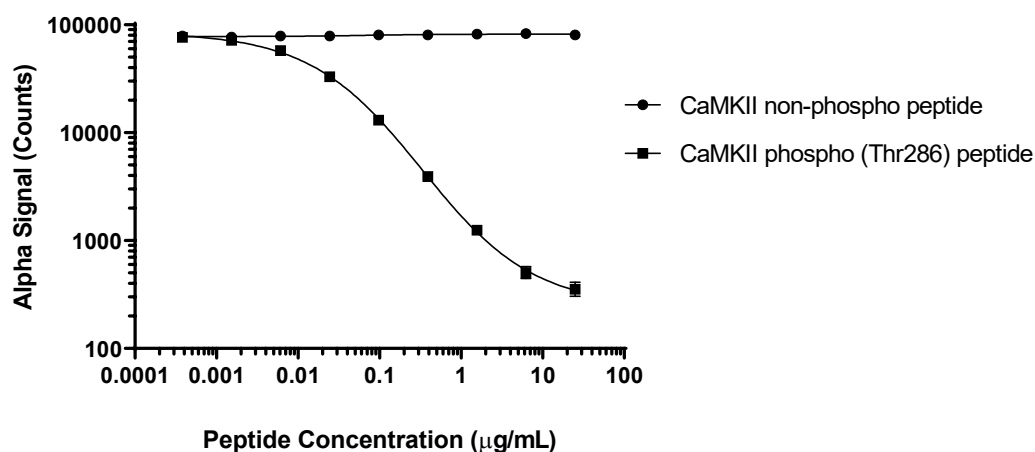
This assay kit contains antibodies which recognize the phospho-Thr286 epitope and a distal epitope on CaMKII. The proteins detected by this kit corresponds to UniProt ID Q9UQM7, Q13554, Q13555, Q13557. CaMKII is also known as calcium/calmodulin dependent protein kinase II. These antibodies recognize CaMKII- α , CaMKII- β , CaMKII- γ and CaMKII- δ of human and mouse origin. Other species should be tested on a case-by-case basis.

Control lysate information:

Positive Control Lysate: Prepared from U-87 MG cells, cultured to confluence in T175 flasks in 10% FBS containing media and lysed with 5 mL of Lysis Buffer.

Representative data

A fixed concentration of SH-SY5Y positive control lysate was spiked with a dose range of phospho (Thr286) and non-phospho CaMKII peptide. Lysates were assayed for Phospho CaMKII (Thr286) in a 384 well assay plate following the SureFire Ultra assay protocol. Equivalent to approximately 12,000 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"