

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ p-PKCθ (Thr538) Assay Kit

**Products: ALSU-PPKC-A500, ALSU-PPKC-A10K, ALSU-PPKC-A50K,
ALSU-PPKC-A-HV**

Kit-Specificity Information

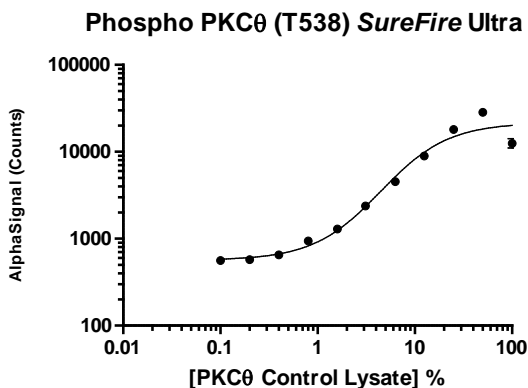
This assay kit contains antibodies which recognize the phospho-Thr538 epitope, and a distal epitope, on PKC-theta (PKCθ). The protein detected by this kit corresponds to Entrez-Gene Id 5588. These antibodies recognize PKCθ of human origin, and predicted to recognise rat and mouse origin protein. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from Jurkat cells, cultured in a T175 flask in 10% FBS containing medium, and lysed at approximately 4×10^6 cells/mL with 1X *Surefire Ultra* Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. Cell lysates, prepared as described above, were diluted with 1X Lysis buffer and assayed.



For more information and detailed protocols, please refer to the generic AlphaLISA *SureFire Ultra* Assay Kits Manual, available from <http://www.perkinelmer.com/category/alpha-surefire-kits>