

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ p-DAP12 (pan Tyr) Assay Kit

ALSU-PDAP12-A500, ALSU-PDAP12-A10K, ALSU-PDAP12-A50K, ALSU-PDAP12-A-HV

Kit Specificity

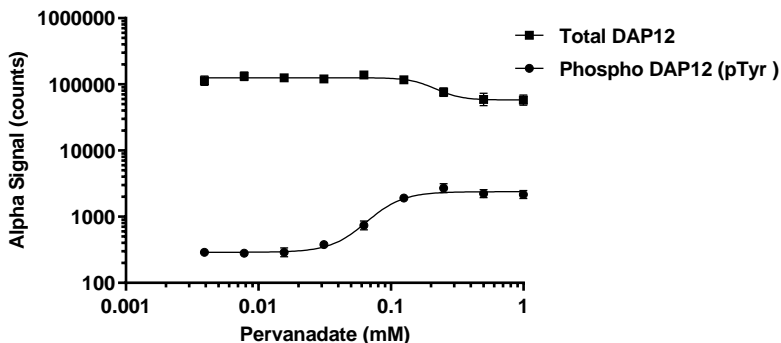
This assay kit contains antibodies which recognize the phospho-pan Tyr, and a distal epitope on DAP12. The protein detected by this kit corresponds to UniProt ID O43914. DAP12 is also known as TYRO protein tyrosine kinase-binding protein and DNAX-activation protein 12. These antibodies recognize DAP12 of human and mouse origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from THP-1 cells cultured to between 0.6-0.8 X10⁶ cells/mL in 10% FBS containing medium. Cells were washed, then treated with 0.25 mM Pervanadate in HBSS for 60 minutes and lysed to 4x10⁶ cells/mL with 5X *SureFire Ultra* Lysis buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. THP1 cells were plated at 600K cells per well in a 96 well plate in HBSS + 0.1% BSA, then treated with Pervanadate at the indicated concentrations for 1 hour at 37°C. Cells were then lysed with 5X Lysis buffer, and assayed for Phospho (pan Tyr) or Total DAP12 in separate wells of an assay plate with respective *SureFire Ultra* kits. Equivalent to approximately 24,000 cells/datapoint.



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>.