

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ DAP12 Total Assay Kit

ALSU-TDAP12-A500, ALSU-TDAP12-A10K, ALSU-TDAP12-A50K, ALSU-TDAP12-A-HV

Kit Specificity

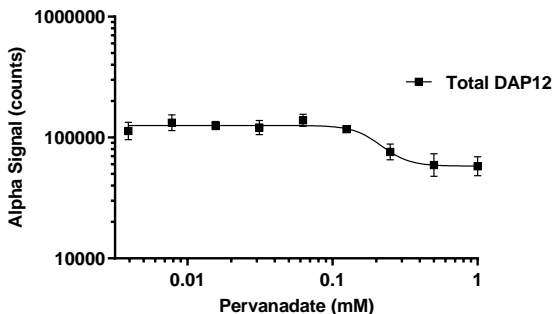
This assay kit contains antibodies which recognize invariant epitopes on DAP12. The protein detected by this kit corresponds to UniProt ID O43914. DAP12 is also known as TYRO protein tyrosine kinase-binding protein and DNAX-activation protein 12. These antibodies recognize DAP12 of human origin. Other 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 and resuspended in HBSS, then 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 Total or Phospho (pan Tyr) 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>.