

AlphaLISA® SureFire® Ultra™

Human DAP12 Total Detection Kit

Product number: ALSU-TDAP12-A500, ALSU-TDAP12-A10K,
ALSU-TDAP12-A50K, ALSU-TDAP12-A-HV



Kit specificity:

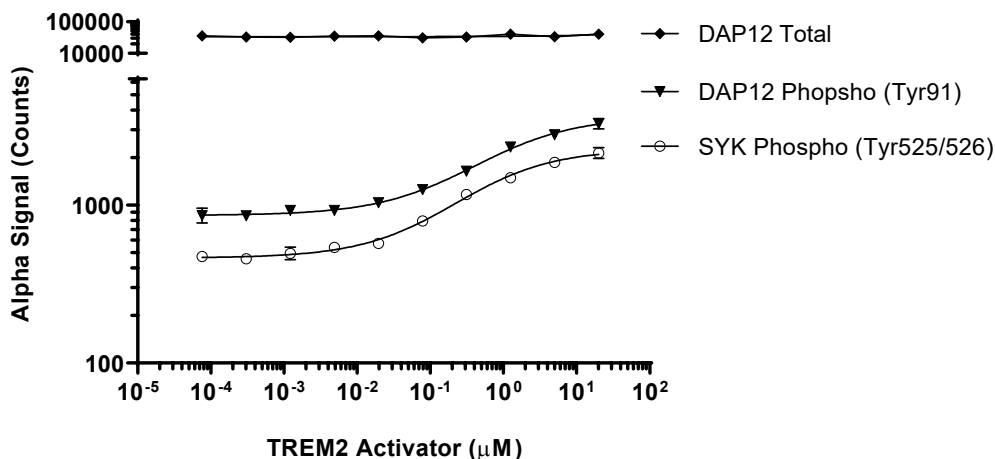
This assay kit contains antibodies which recognize distinct epitopes on DAP12. The protein detected by this kit corresponds to UniProt ID O43914. DAP12 is also known as TYRO protein tyrosine kinase-binding protein and DNAX-activation protein 12. These antibodies recognize DAP12 of human origin. Other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

Control lysate information:

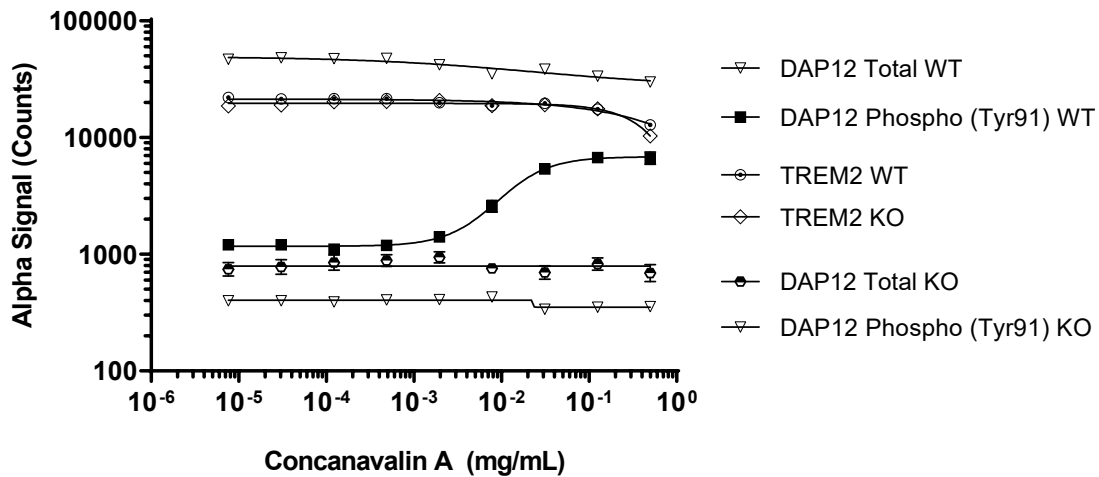
Positive Control Lysate: Prepared from THP 1 cells cultured in 10% FBS containing media. Cells were harvested and washed with HBSS + 0.1% BSA, adjusted to 4 x10⁶ cells/mL and then treated with 0.5 mM Pervanadate for 30 minutes. Cells were then lysed with the addition of 5X Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. THP1 cells were prepared at 600K/mL in 10% FBS containing media and treated with 35 ng/mL TGFβ for 20 hours. Cells were then washed with HBSS + 0.1% BSA and seeded at 400K cells/well in a 96 well plate. Cells were treated with a TREM2 Activator at the indicated concentrations for 10 minutes and lysed with 5X Lysis Buffer. Assayed separately for Phospho (Tyr91) and Total DAP12, and Phospho SYK (Tyr525/526) using respective *SureFire Ultra* kits. Equivalent to approximately 16,000 cells or 1,600 cells/datapoint (DAP12 Total).



Data obtained with a 2-plate, 2-incubation protocol. THP1 Wild Type cells (WT) and THP1-DAP12 Knockout cells (KO) were harvested at 800K/mL and washed in HBSS + 0.1% BSA. Cells were seeded at 400K cells/well in a 96 well plate and incubated with Concanavalin A at the indicated concentrations for 10 minutes. Cells were spun down, washed in HBSS and lysed in Lysis Buffer. Assayed separately for Phospho (Tyr91), Total DAP12 and Total TREM2 using respective *SureFire Ultra* kits. Equivalent to approximately 20,000 cells/datapoint (Phospho DAP12 Tyr91) or 2,000 cells/datapoint (Total DAP12 and Total TREM2).



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"

