

AlphaLISA™ SureFire® Ultra™

Human p-SIRPα (pan Tyr) Detection Kit

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



Kit specificity:

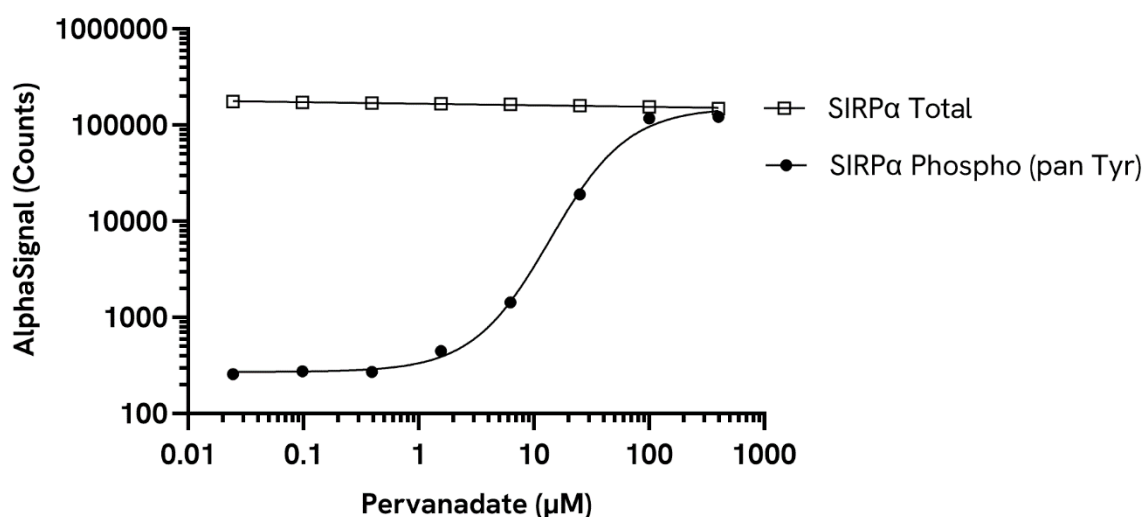
This assay kit contains antibodies which recognize phospho-Tyr and a distal epitope on SIRPα. The protein detected by this kit corresponds to UniProt ID P78324. SIRPα is also known as SHPS1, MYD1 and PTPNS1. These antibodies recognize SIRPα of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

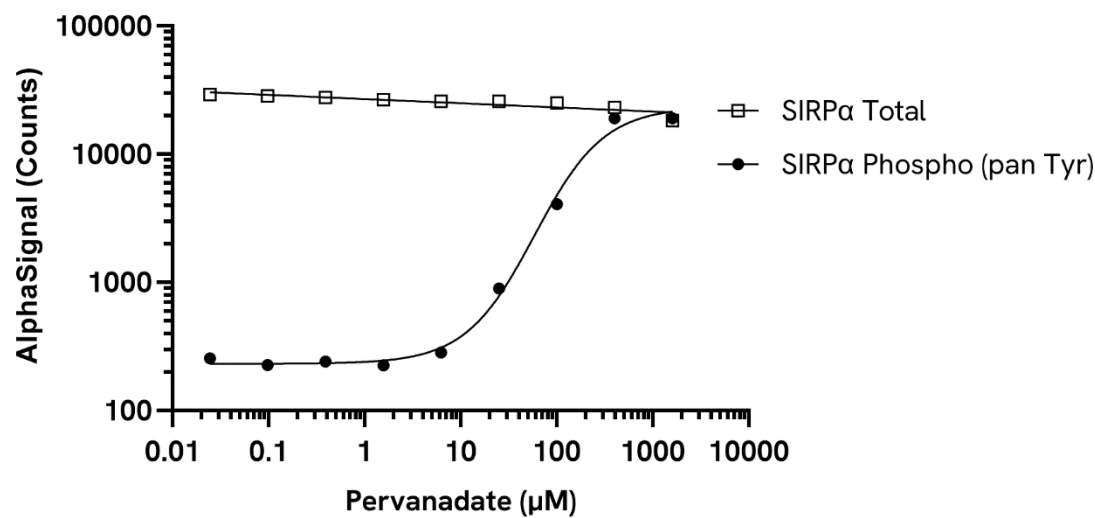
Positive Control Lysate: Prepared from U937 cells cultured in 10% FBS containing medium. Cells were harvested, washed with HBSS, adjusted to 10⁶ cells/mL and treated with 500 μM Pervanadate for 15 minutes. Cells were lysed with the addition of 5X Lysis Buffer.

Representative data:

Data obtained with a 2-plate, 2-incubation protocol. U937 cells were washed in HBSS + 0.1% BSA, seeded at 50K cells/well in a 96-well plate and treated with Pervanadate at the indicated concentrations for 15 minutes. Cells were lysed with 5X Lysis Buffer and assayed separately for Phospho (pan Tyr) and Total SIRPα using respective SureFire Ultra kits. Equivalent to approximately 2,000 cells/datapoint.



Data obtained with a 2-plate, 2-incubation protocol. THP-1 cells were washed in HBSS + 0.1% BSA, seeded at 50K cells/well in a 96-well plate and treated with Pervanadate at the indicated concentrations for 15 minutes. Cells were lysed with 5X Lysis Buffer and assayed separately for Phospho (pan Tyr) and Total SIRPα using respective *SureFire Ultra* kits. Equivalent to approximately 2,000 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"

