

AlphaLISA® SureFire® Ultra™

Human p-PD-1 (pan Tyr) Detection Kit

Product number: ALSU-PPD1-A500, ALSU-PPD1-A10K,
 ALSU-PPD1-A50K, ALSU-PPD1-A-HV



Kit specificity:

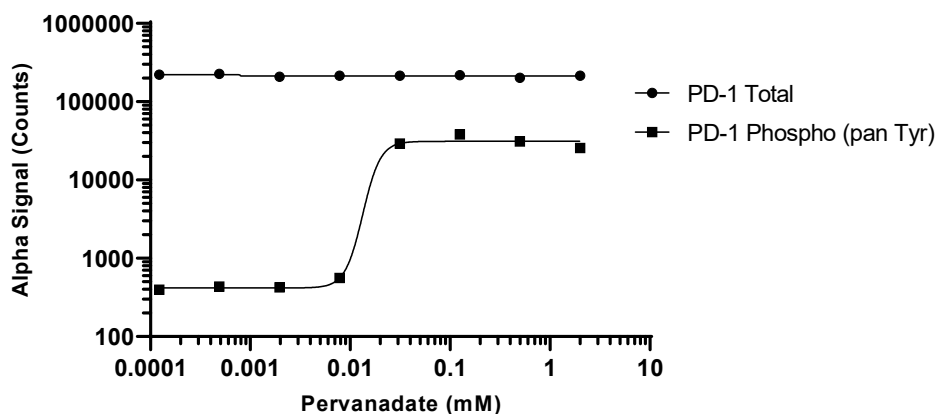
This assay kit contains antibodies which recognize phospho-tyrosine epitopes and an invariant epitope on programmed cell death 1 protein (PD-1). The protein detected by this kit corresponds to UniProt ID Q15116. PD-1 is also known as PDCD1. These antibodies recognize PD-1 of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

A control lysate is not available for this target. However, PD-1 CHO cells, in which the PD-1 protein is transfected, are available from Abeomics (Product code: 14-500ACL).

Representative data

Data obtained with a 2-plate, 2-incubation protocol. PD-1 CHO cells were plated at 40K cells/well in a 96 well plate and incubated overnight. Cells were then treated with Pervanadate at the indicated concentrations for 30 minutes. Cells were lysed with Lysis Buffer and assayed for Phospho-tyrosine and Total PD-1 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"