

# Technical Data Sheet

## AlphaLISA® SureFire® Ultra™ p-PD-1 (pan Tyr) Assay Kit

ALSU-PPD1-A500, ALSU-PPD1-A10K, ALSU-PPD1-A50K, ALSU-PPD1-A-HV

### Kit-Specificity

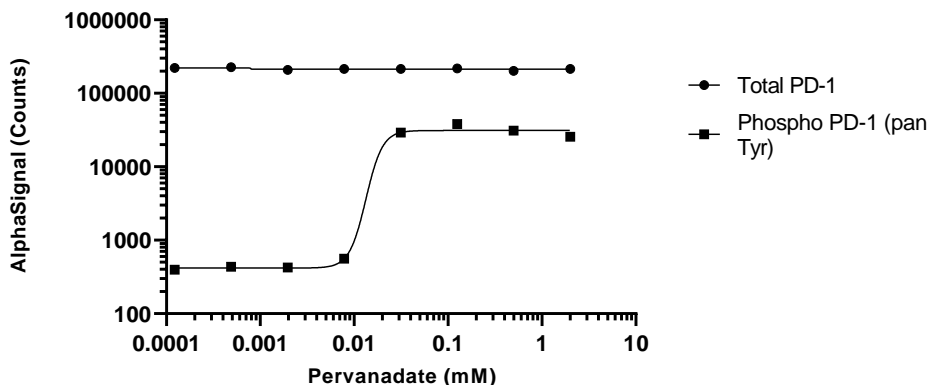
This assay kit contains antibodies which recognize phosphotyrosine epitopes and an invariant epitope on programmed cell death 1 protein (PD-1). The protein detected by this kit corresponds to UniProt ID Q15116. PD-1 is also known as PDCD1. These antibodies recognize PD-1 of human origin. Other species 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 per 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 1x Lysis Buffer and assayed for Total or Phospho PD-1 (pan-Tyr) using respective *SureFire Ultra* kits. Equivalent to approximately 2000 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>