

Technical Data Sheet

Caution: Research use only.

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

**Products: ALSU-PPD1-A500, ALSU-PPD1-A10K, ALSU-PPD1-A50K,
ALSU-PPD1-A-HV**

Kit-Specificity Information

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 GenBank Accession NG_012110.1. PD-1 is also known as PDCD1 or CD279. These antibodies recognize PD-1 of human origin. Other species should be tested on a case-by-case basis.

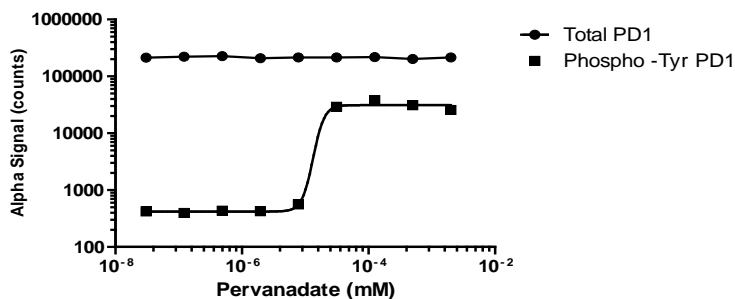
Control Lysate Information

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 seeded at 20K cells per well, grown overnight, and then treated with the indicated concentrations of Pervanadate for 30min, and lysed with *SureFire Ultra* 1X Lysis buffer. Samples were assayed for phospho PD-1 (pan-Tyr) and Total PD-1, using respective *SureFire Ultra* kits.

Pervanadate Dose Response curves in PD1 CHO cells PD-1 Ultra assays



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>