

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ p-SHP-1 (Tyr536) Assay Kit

ALSU-PSHP1-A500, ALSU-PSHP1-A10K, ALSU-PSHP1-A50K, ALSU-PSHP1-A-HV

Kit-Specificity

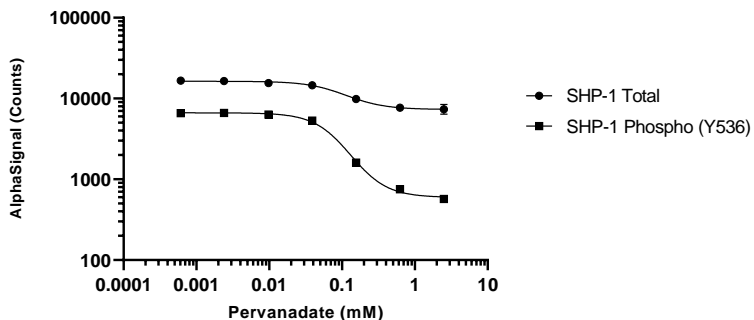
This assay kit contains antibodies that specifically recognize the phospho-Tyr536-epitope and a distal epitope on SHP-1. The protein detected by this kit corresponds to UniProt ID P29350. SHP-1 is also known as Tyrosine-protein phosphatase non-receptor type 6, hematopoietic cell protein-tyrosine phosphatase, PTP-1C and SH-PTP1. These antibodies recognize SHP-1 of human origin. Rat and mouse are not detected. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from Jurkat cells cultured in T175 flasks to 1×10^6 cells/mL in medium containing 10% FBS and lysed at 3.2×10^6 cells/mL with 1X *SureFire Ultra* Lysis Buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. Jurkat cells were washed and resuspended in HBSS at a density of 2×10^6 cells/mL. Cells were treated with Pervanadate at the indicated concentrations for 30 minutes and lysed with 5x Lysis Buffer at a final cell density of 0.8×10^6 cells/mL. Cell lysates were assayed separately for Total or Phospho (Tyr536) SHP-1 using respective *SureFire Ultra* kits. Equivalent to approximately 8000 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>