

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ p-SHP-1 (Tyr564) Assay Kit

**Products: ALSU-PSHP1-B500, ALSU-PSHP1-B10K, ALSU-PSHP1-B50K,
ALSU-PSHP1-B-HV**

Kit-Specificity Information

This assay kit contains antibodies that specifically recognize the phospho-Tyr564-epitope, or an invariant epitope on SHP-1. The human protein detected by this kit corresponds to Entrez Gene ID 5777 (human).

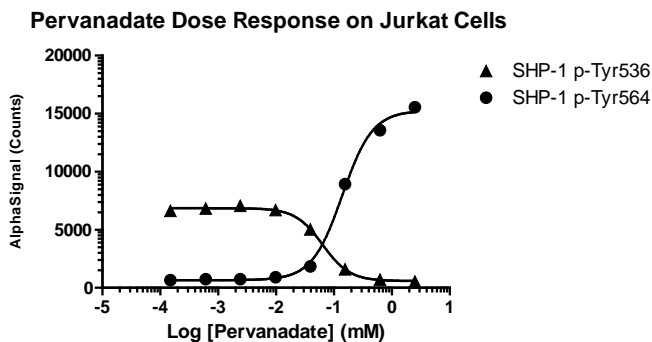
These antibodies recognize SHP-1 of human origin. Rat and mouse are not detected. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from Jurkat cells cultured in a T175 flask to 10^6 cells/mL in medium containing 10%FBS, treated with 2.5mM Pervanadate for 30 minutes, 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 (0.2×10^6 cells/well) stimulated for 30min with Pervanadate. Cells were lysed with 5X Lysis buffer in culture medium. Data shows phosphorylation of both SHP-1 Tyr536 and SHP-1 Tyr564 (measured with respective kits).



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>