

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ p-SLP-76 (Tyr113) Assay Kit

ALSU-PSLP-B500, ALSU-PSLP-B10K, ALSU-PSLP-B50K, ALSU-PSLP-B-HV

Kit Specificity

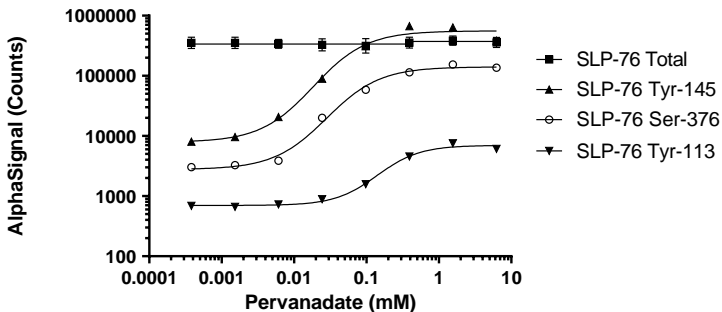
This assay kit contains antibodies which recognize the phospho-Tyr113 epitope, and a distal epitope, on SLP-76. The protein detected by this kit corresponds to UniProt ID 13094. SLP-76 is also known as lymphocyte cytosolic protein 2. These antibodies recognize SLP-76 of human origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from Jurkat cells, cultured in T175 flask to 10^6 cells/mL in 10% FBS containing medium, then treated with 11 mM H_2O_2 and 50 nM Calyculin A for 10 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 harvested from a flask at 500 K/mL were seeded at 4×10^6 cells/mL in HBSS + 0.1% BSA in a 96 well plate. Cells were treated with a dose range of pervanadate for 30 mins before lysing with 5X *SureFire Ultra* Lysis buffer. Lysates were assayed for Phospho (Tyr-113, Tyr-145 & Ser-376) and Total SLP-76 in separate wells of an assay plate using respective *SureFire Ultra* kits. Equivalent to approximately 16,000 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>.