

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ SLP-76 Total Assay Kit

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

Kit Specificity

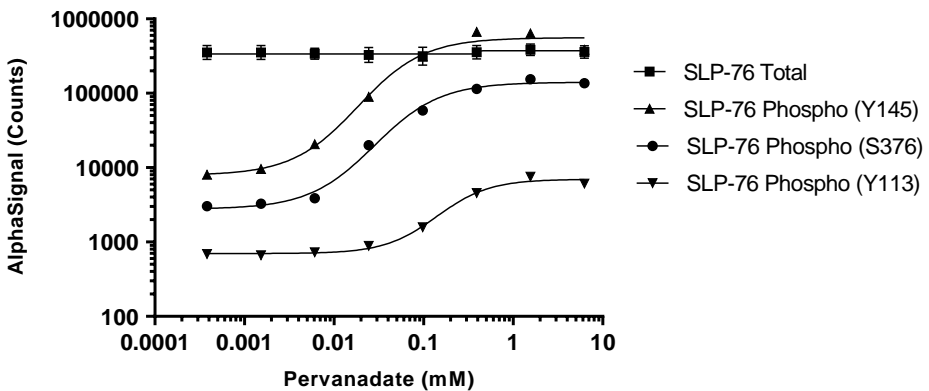
This assay kit contains antibodies which recognize invariant epitopes on SLP-76. The protein detected by this kit corresponds to UniProt ID Q13094. SLP-76 is also known as lymphocyte cytosolic protein 2. These antibodies recognize SLP-76 of human origin. Other species should be tested on a case-by-case 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 were harvested, washed and resuspended in HBSS + 0.1% BSA at a density of 4×10^6 cells/mL. Cells were then treated with pervanadate at the indicated concentrations for 30 minutes. Cells were lysed with 5x Lysis Buffer and assayed separately for Total or Phospho SLP-76 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>