

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ p-SYK (Tyr352) Assay Kit

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

Kit-Specificity

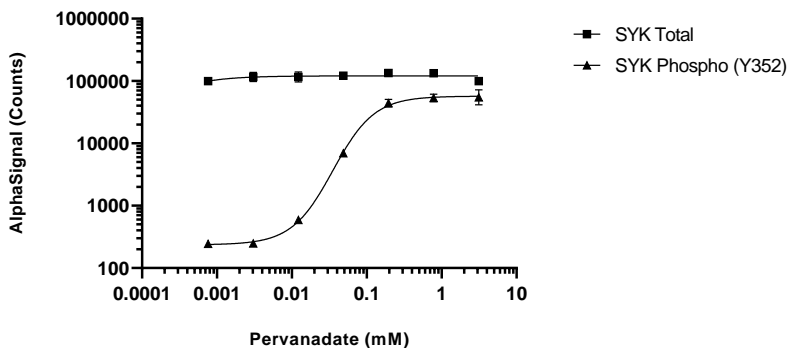
This assay kit contains antibodies which recognize the phospho-Y352 epitope, and distal epitope on Tyrosine-protein kinase SYK. The protein detected by this kit corresponds to UniProt ID P43405. SYK is also known as spleen tyrosine kinase and p72-SYK. These antibodies recognize SYK of human origin. It is predicted to react also with mouse and rat. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from U937 cells, cultured to approximately 1×10^6 cells/mL in T175 flasks in 10% FBS containing medium, then treated with 0.5 mM Pervanadate for 30 minutes and lysed at 1×10^6 cells/mL with 1X *SureFire Ultra* Lysis Buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. U937 cells were harvested, washed in HBSS and resuspended at a density of 2×10^6 cells/mL, then treated with Pervanadate at the indicated concentrations for 20 minutes. Cells were lysed with 1x Lysis Buffer at a final cell density of 0.8×10^6 cells/mL and assayed for Total or Phospho (Y525/526) SYK 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>