

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ p-SYK (Tyr525/526) Assay Kit

**Products: ALSU-PSYK-A500, ALSU-PSYK-A10K, ALSU-PSYK-A50K,
ALSU-PSYK-A-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Tyr525/526 epitope, and distal invariant epitopes, on SYK. The protein detected by this kit corresponds to Entrez-Gene ID 6850.

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

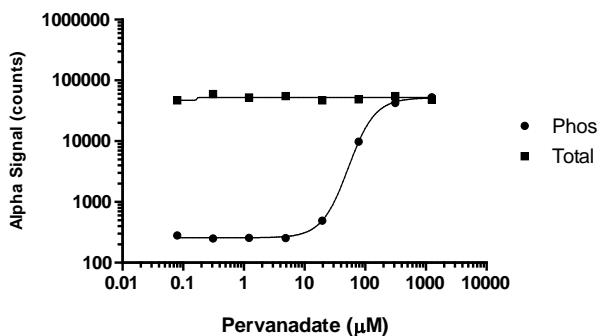
Positive Control Lysate: Prepared from U937 cells, cultured to 10^6 cells/ml in T175 flasks in 10% FBS containing medium, then treated with 1% Pervanadate for 30min and lysed to 10^6 cells/mL with 1X *SureFire Ultra* Lysis buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol.

Phospho and Total SYK were then measured in separate assay wells using *SureFire Ultra* assay kits.

Pervanadate stimulation of p-SYK in U937 cells



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>