

# Technical Data Sheet

Caution: Research use only.

## AlphaLISA® SureFire® Ultra™ SYK Total Assay Kit

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

### **Kit-Specificity Information**

This assay kit contains antibodies which recognize two 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**

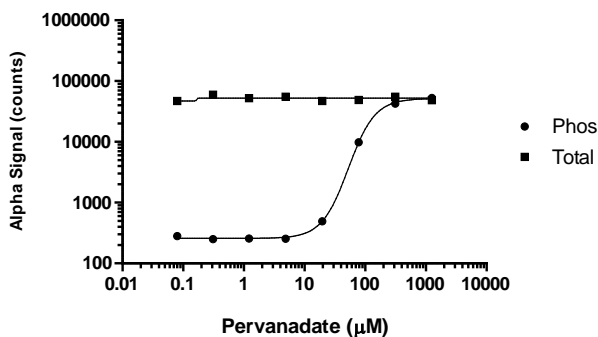
Positive Control Lysate: Prepared from U937 cells, cultured to  $10^6$  cells/ml in T175 flasks in 10% FBS containing medium, then treated with 1% Pervanadate for 30min and lysed to  $10^6$  cells/mL with 1X *SureFire Ultra* Lysis buffer.

### **Representative data**

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

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

#### **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>