

# Technical Data Sheet

Caution: Research use only.

## Alpha SureFire® Ultra™ Multiplex p-SYK (Tyr525/526) + Total SYK Assay Kit

**Products: MPSU-PTSYPK-K500, MPSU-PTSYPK-K10K, MPSU-PTSYPK-K50K,  
MPSU-PTSYPK-K-HV**

### **Kit-Specificity Information**

This assay kit contains antibodies which recognize the phospho-Tyr525/526 epitope, and a distal epitope, on SYK for phospho SYK analysis, plus antibodies to measure the total levels of SYK.

These antibodies recognize SYK of human origin. They are predicted to react also with mouse and rat. Other species should be tested on a case-by-case basis.

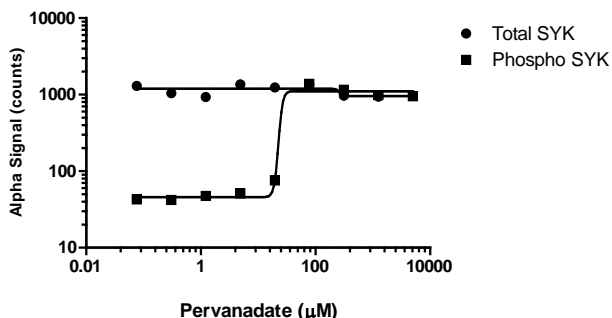
### **Control Lysate Information**

Positive Control Lysate: Prepared from U937 cells, cultured to  $10^6$  cells/ml in T175 flasks in 10% FBS containing medium, then treated with 1mM 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. U937 cells were plated at 200K cells per well in a 96 well plate, and treated with Pervanadate for 30 minutes at the indicated concentrations. Cells were lysed in medium with 5X Lysis buffer and assayed for both phospho and total SYK as detailed in the MPSU manual.

#### **Pervanadate Dose Response in U937 cells - SYK Multiplex**



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>