

# Technical Data Sheet

Caution: Research use only.

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

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

### **Kit-Specificity Information**

This assay kit contains antibodies which recognize two invariant epitopes on SLP-76. The protein detected by this kit corresponds to GenBank Accession U20158.1. These antibodies recognize SLP-76 of human origin. Other species should be tested on a case-by-case basis.

### **Control Lysate Information**

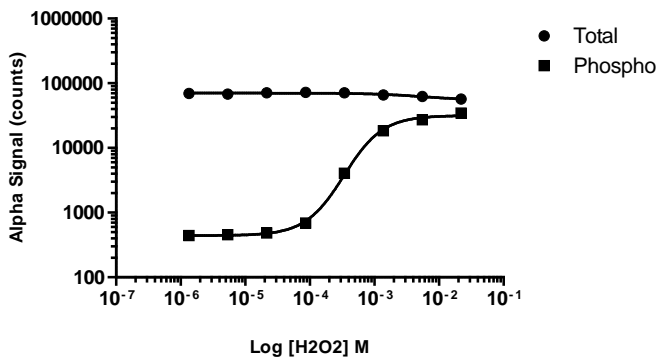
Positive Control Lysate: Prepared from Jurkat cells, cultured in T175 flask to  $10^6$  cells/mL in 10% FBS containing medium, then treated with 11 mM  $H_2O_2$  and 50  $\mu$ M Calyculin for 10 mins and lysed at  $3.2 \times 10^6$  cells/mL with 1X *SureFire Ultra* Lysis buffer.

### **Representative data**

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

Lysates were assayed for p-SLP-76 and total SLP-76 in separate wells using the respective *SureFire Ultra* kits.

#### **Total-Phospho SLP-76 - Jurkat 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>