

# Technical Data Sheet

Caution: Research use only.

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

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

### **Kit-Specificity Information**

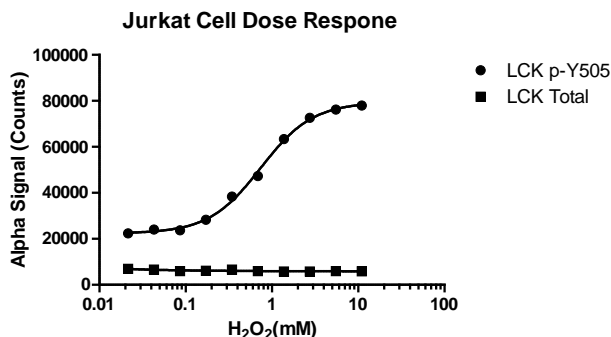
This assay kit contains antibodies which recognize two invariant epitopes on Lck. The protein detected by this kit corresponds to GenBank Accession NG\_023387.1. These antibodies recognize Lck of human, mouse and rat origin. Other species should be tested on a case-by-case basis.

### **Control Lysate Information**

Positive Control Lysate: Prepared from Jurkat cells, cultured in a T175 flask to approximately  $3 \times 10^6$  cells/mL in 10% FBS containing medium, then treated with 11mM Peroxide plus  $50 \mu\text{M}$  Calyculin A for 10min and lysed with 5X *SureFire Ultra* Lysis buffer.

### **Representative data**

Data obtained with a 2-plate, 2-incubation protocol. Jurkat cells were plated at 100K cells per well, and treated with Peroxide at the indicated concentrations for 10 minutes, then lysed in medium with 5X Lysis buffer. Levels of Lck (Tyr505) phosphorylation, and Total Lck were assayed with respective *SureFire Ultra* assay kits.



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>