

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ p-Lck (Tyr505) Assay Kit

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

Kit-Specificity Information

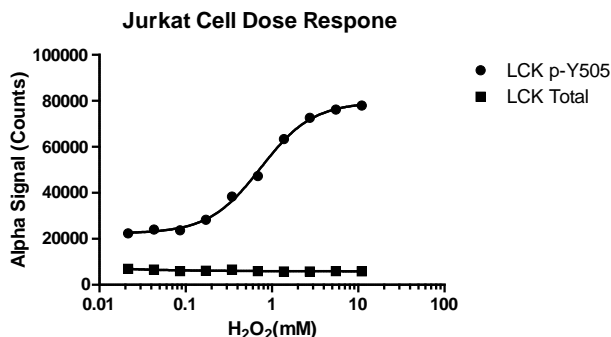
This assay kit contains antibodies which recognize the phospho-Tyr505 epitope, and a distal epitope, on Lck. The protein detected by this kit corresponds to GenBank Accession NG_023387.1. These antibodies recognize Lck of human origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from Jurkat cells, cultured in a T175 flask to approximately 3×10^6 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>