

# Technical Data Sheet

## AlphaLISA® SureFire® Ultra™ PLCγ1 Total Assay Kit

ALSU-TPLCG1-A500, ALSU-TPLCG1-A10K, ALSU-TPLCG1-A50K, ALSU-TPLCG1-A-HV

### Kit Specificity

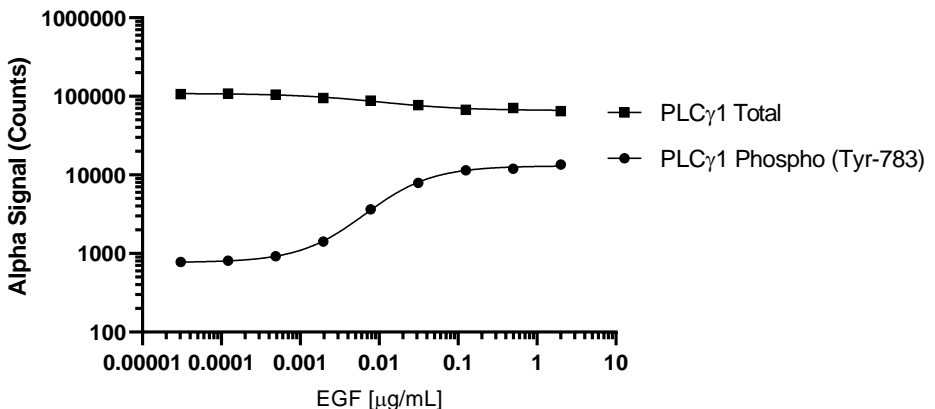
This assay kit contains antibodies which recognize invariant epitopes on PLCγ1. The protein detected by this kit corresponds to UniProt ID P19174. These antibodies recognize PLCγ1 of human and mouse origin. Other species should be tested on a case-by-case basis.

### Control Lysate Information

Positive Control Lysate: Prepared from A431 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 250 ng/mL EGF for 10 minutes and lysed in 4 mL of 1X *SureFire Ultra* Lysis buffer.

### Representative Data

Data obtained with a 2-plate, 2-incubation protocol. A431 cells were seeded overnight at 40K cells/well in a 96 well plate in DMEM + 10% FBS. Cells were treated with EGF at the indicated concentrations for 10 minutes at 37°C. Cells were then lysed with 1X *SureFire Ultra* Lysis buffer and assayed for Phospho (Tyr783) and Total PLCγ1 in separate wells of an assay plate using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint.



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>.