

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ p- PLCγ1 (Ser1248) Assay Kit

ALSU-PPLCG1-B500, ALSU-PPLCG1-B10K, ALSU-PPLCG1-B50K, ALSU-PPLCG1-B-HV

Kit Specificity

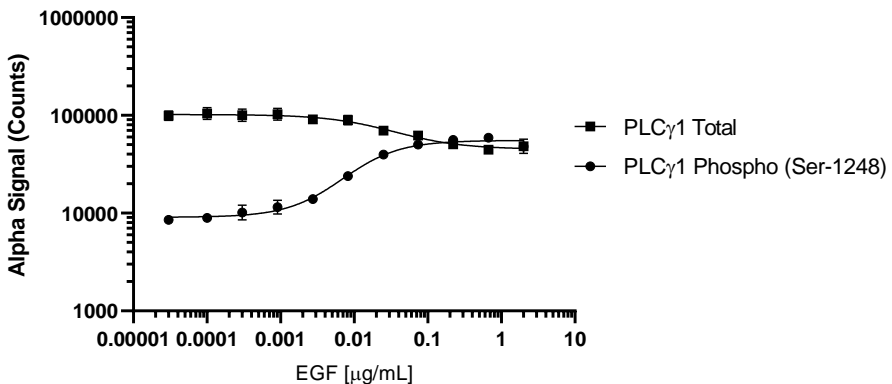
This assay kit contains antibodies which recognize phospho Ser-1248 and a distal epitope on PLCγ1. The protein detected by this kit corresponds to UniProt ID P19174. These antibodies recognize PLCγ1 of human and mouse origin. Other species should be tested on a case-by-case basis.

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 (Ser1248) 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>