

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ p-EGF Receptor (Tyr1068) Assay Kit

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

Kit-Specificity

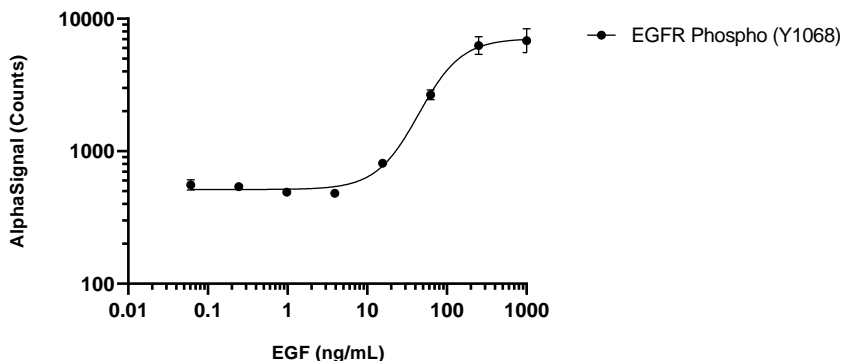
This assay kit contains antibodies which recognize the phospho-Tyr1068 epitope, and a distal epitope, on epidermal growth factor receptor (EGFR). The protein detected by this kit corresponds to UniProt ID P00533. EGFR is also known as Proto-oncogene c-ErbB-1 and Receptor tyrosine-protein kinase erbB-1. These antibodies recognize EGFR 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 2 µg/mL EGF for 10 minutes and lysed in 4 mL of 1X Lysis Buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. A431 cells were plated at 40K cells per well in a 96 well plate and incubated overnight. Cells were serum starved for 2 hours prior to treatment with EGF at the indicated concentrations for 10 minutes. Cells were lysed with 1x Lysis Buffer and assayed for Total or Phospho (Tyr1068) EGFR with respective SureFire Ultra kits. Equivalent to approximately 200 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>