

# Technical Data Sheet

## AlphaLISA® SureFire® Ultra™ EGF Receptor Total Assay Kit

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

### Kit-Specificity

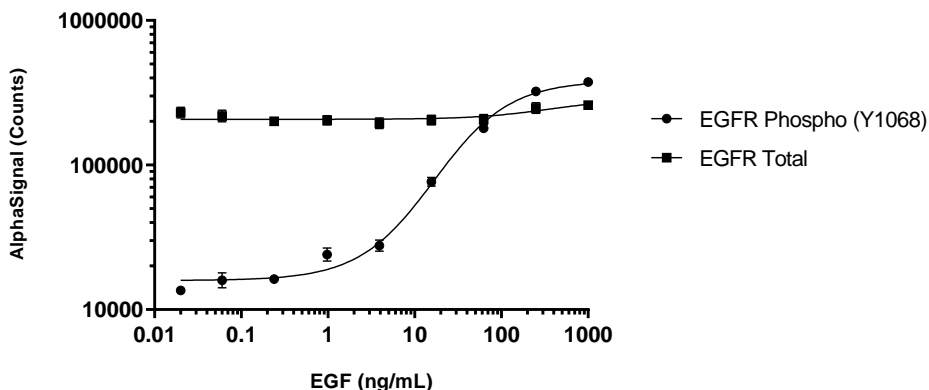
This assay kit contains antibodies which recognize invariant epitopes on EGF Receptor. The protein detected by this kit corresponds to UniProt ID P00533. EGFR is also known as Proto-oncogene c-ErbB-1 and Receptor tyrosine-protein kinase erbB-1. These antibodies recognize EGF Receptor of human and mouse 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 10min and lysed in 4mL of 1X *SureFire Ultra* 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 400 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>