

# Technical Data Sheet

Caution: Research use only.

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

**Products: ALSU-PEGFR-B500, ALSU-PEGFR-B10K, ALSU-PEGFR-B50K,  
ALSU-PEGFR-B-HV**

### **Kit-Specificity Information**

This assay kit contains antibodies which recognize the phospho-Tyr1068 epitope, and a distal epitope, on EGF Receptor. The protein detected by this kit corresponds to GenBank Accession NP\_005219. EGFR is also known as ERBB, HER1, mENA, ERBB1 and PIG61. These antibodies recognize EGF Receptor of human and mouse origin. Other species should be tested on a case-by-case basis.

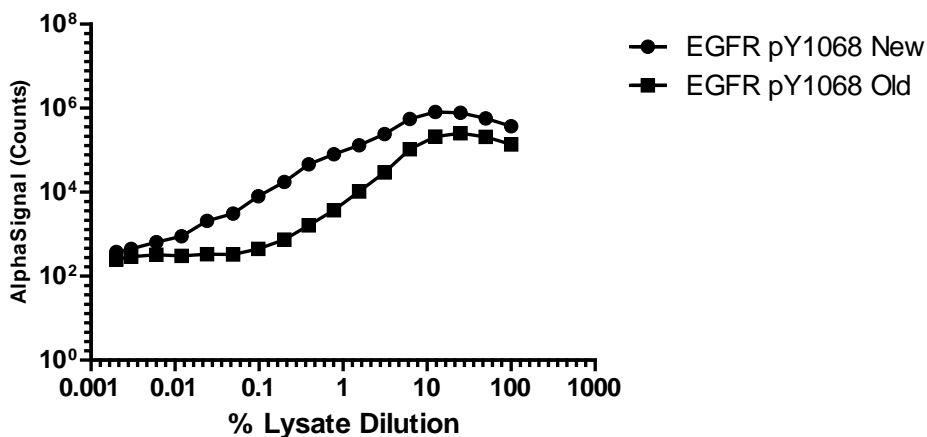
### **Control Lysate Information**

Positive Control Lysate: Prepared from A431 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 2µg/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. Control Lysate was diluted as indicated in 1X Lysis Buffer, and assayed according to the protocol.

#### **Lysate Dilution of EGFR pY1068 New vs Old**



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>