

Technical Data Sheet

Caution: Research use only.

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

**Products: ALSU-PEGFR-A500, ALSU-PEGFR -A10K, ALSU-PEGFR-A50K,
ALSU- PEGFR-A-HV**

Kit-Specificity Information

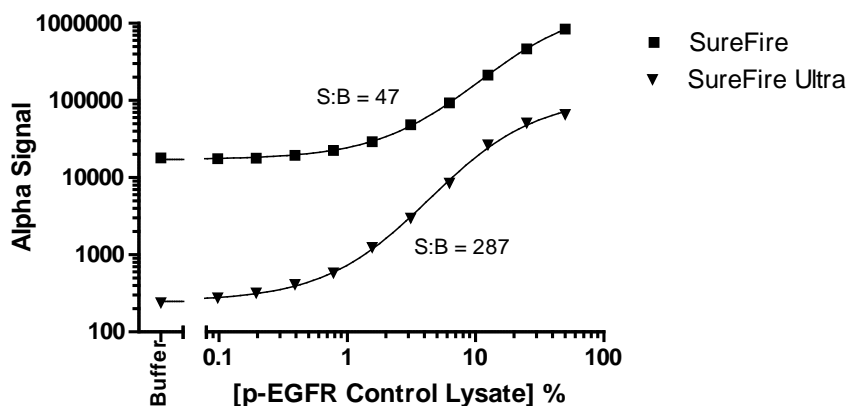
This assay kit contains antibodies which recognize the phospho-Tyr1068 epitope, and a distal epitope, on epidermal growth factor receptor (EGFR). The protein detected by this kit corresponds to GenBank Accession NP_005219. EGFR is also known as ERBB, HER1, mENA, ERBB1 and PIG61. These antibodies recognize 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 2µg/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. Data shows a comparison of AlphaScreen SureFire and AlphaLISA SureFire Ultra assay performance. Control lysate was diluted in the Lysis buffer specific for the assay type.



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>