

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex p-EGF Receptor (Tyr1068) + Total EGF Receptor Assay Kit

**Products: MPSU-PTEGFR-K500, MPSU-PTEGFR-K10K, MPSU-PTEGFR-K50K,
MPSU-PTEGFR-K-HV**

Kit-Specificity Information

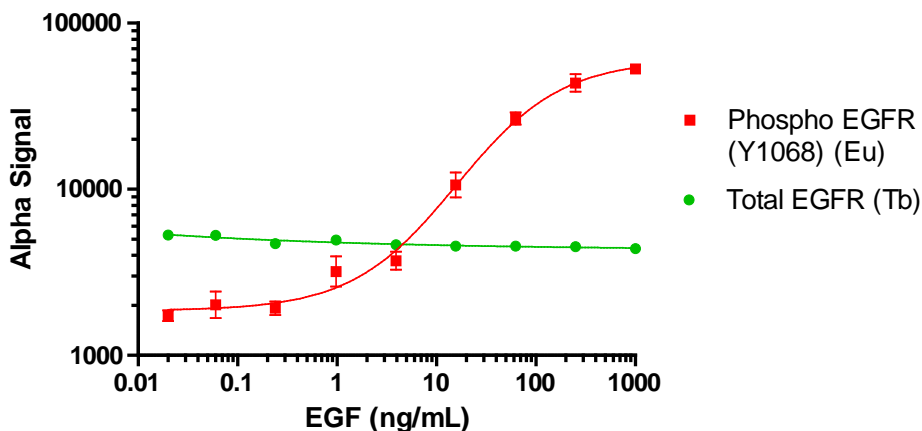
This assay kit contains antibodies which recognize the phospho-Tyr1068 epitope, and a distal epitope, on EGF Receptor (EGFR) for phospho EGFR analysis, plus antibodies to measure the total levels of EGFR. These antibodies recognize EGF Receptor of human and mouse origin. Other species should be tested 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 μ g/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 200K cells per mL, incubated overnight, starved for 2 hours, then stimulated with EGF for 10 minutes at the indicated concentrations. Cells were lysed with 100 μ L of 1X Lysis buffer and assayed for both phospho and total EGFR as detailed in the MPSU manual.



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>