

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ EGF Receptor Total Assay Kit

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

Kit-Specificity Information

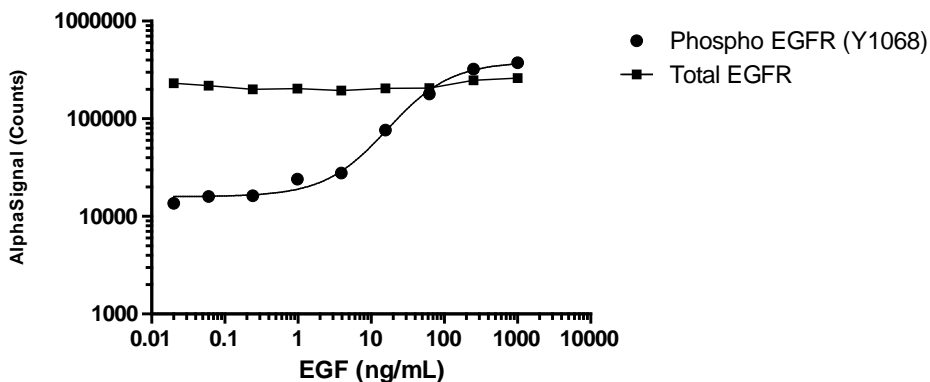
This assay kit contains antibodies which recognize invariant epitopes on EGF Receptor. These antibodies recognize EGF Receptor of human and mouse origin. Other species should be tested on a case-by-case basis.

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. 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 100µL of 1X Lysis buffer and assayed.



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>