

AlphaLISA® SureFire® Ultra™

Human EGF Receptor Total Detection Kit

Product number: ALSU-TEGFR-A500, ALSU-TEGFR-A10K,
ALSU-TEGFR-A50K, ALSU-TEGFR-A-HV



Kit specificity:

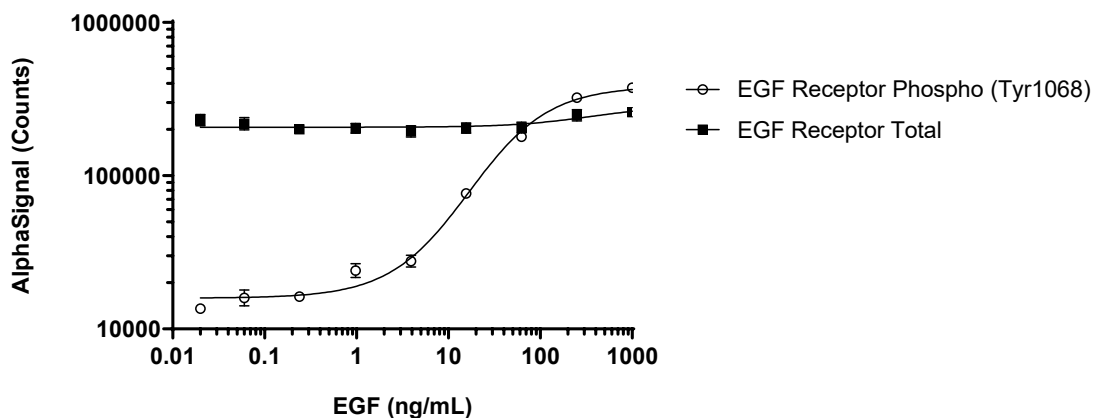
This assay kit contains antibodies which recognize distinct epitopes on EGF Receptor. The protein detected by this kit corresponds to UniProt ID P00533. EGF Receptor is also known as epidermal growth factor receptor, c-ErbB1 and HER1. These antibodies recognize EGFR of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

Positive Control Lysate: Prepared from A-431 cells, cultured to confluence in T175 flasks in 10% FBS containing media, treated with 2 µg/mL EGF for 10 minutes and lysed with 8 mL of Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. A-431 cells were seeded at 40K cells per well in a 96 well plate and incubated overnight. Cells were serum starved for 2 hours and then treated with EGF at the indicated concentrations for 10 minutes. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Tyr1068) or Total EGF Receptor using respective *SureFire Ultra* kits. Equivalent to approximately 400 cells/datapoint.



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF"