

AlphaLISA® SureFire® Ultra™

# Human EGF Receptor L858R Total Detection Kit

**Product number:** ALSU-TEGFR-B500, ALSU-TEGFR-B10K,  
ALSU-TEGFR-B50K, ALSU-TEGFR-B-HV



**Kit specificity:**

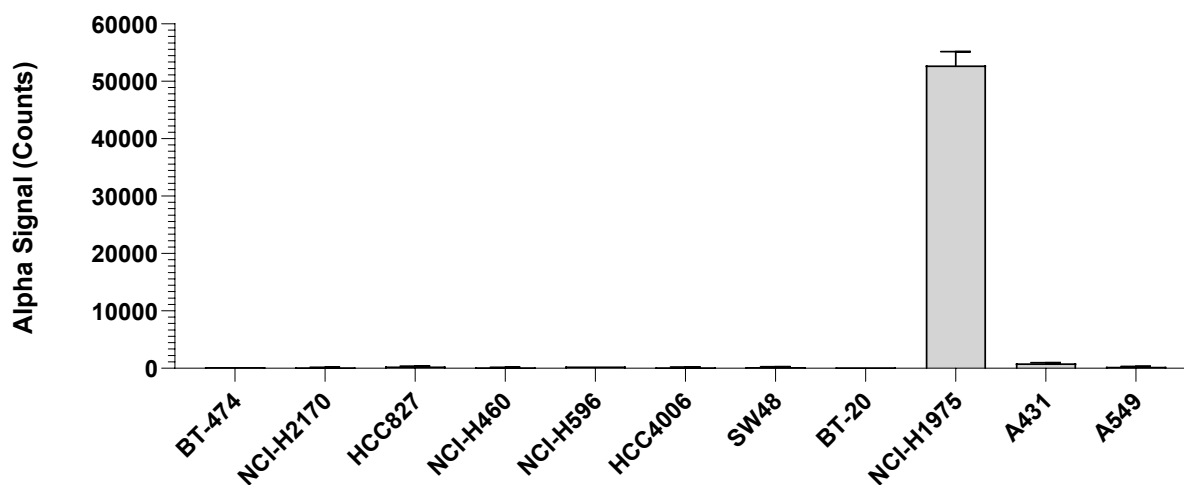
This assay kit contains antibodies which recognize the L858R mutation and a distinct epitope on EGF Receptor. The protein detected by this kit corresponds to UniProt ID P00533. EGF Receptor is also known as EGFR, ERBB1 and HER1. These antibodies recognize EGF Receptor L858R of human origin. Other species should be tested on a case-by-case basis.

**Control lysate information:**

Positive Control Lysate: Prepared from NCI-H1975 cells, cultured to confluence in T175 flasks in 10% FBS containing media and lysed with 4 mL of Lysis Buffer.

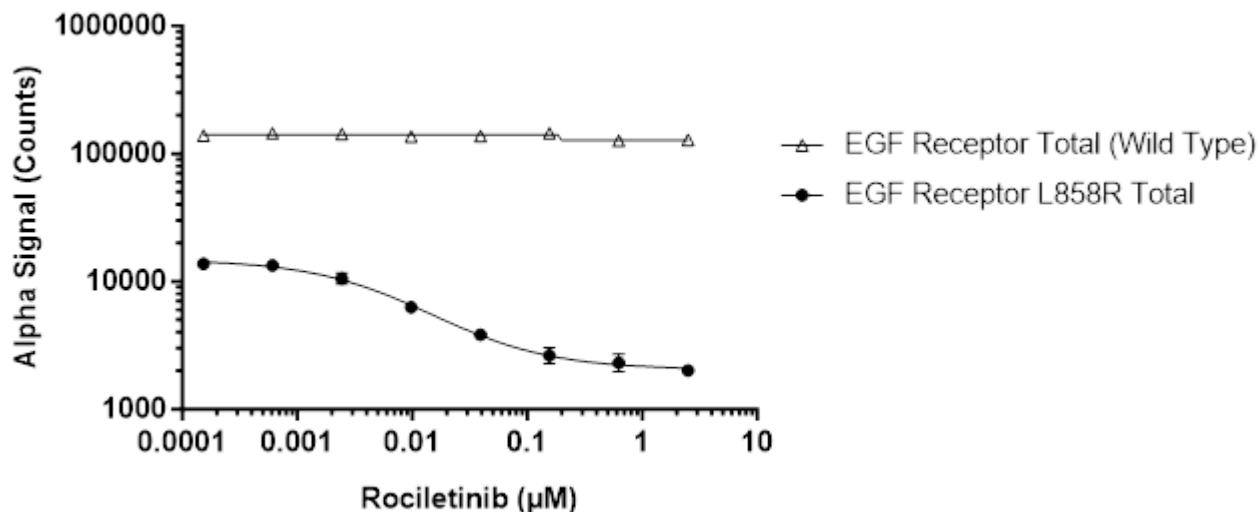
**Representative data**

Data obtained for EGF Receptor L858R Total assay from screening various lysates prepared from cell lines that carry hotspot mutations or gene copy number amplification within the EGFR or ERBB2 gene, in addition to cell lines expressing wild type EGFR (A431, A459). Approximate number of cells/datapoint is indicated for the various cell lines.



Approx. 20,000 cells/datapoint

Data obtained with a 2-plate, 2-incubation protocol. NCI-H1975 cells were seeded at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with Rociletinib (L858R mutant-selective EGFR inhibitor) at the indicated concentrations for 24 hours. Cells were lysed with Lysis Buffer and assayed separately for Total EGFR Receptor L858R and EGFR Receptor (Wild Type) using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 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"

