

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ ErbB2 Total Assay Kit

ALSU-TEB2-A500, ALSU-TEB2-A10K, ALSU-TEB2-A50K, ALSU-TEB2-A-HV

Kit-Specificity

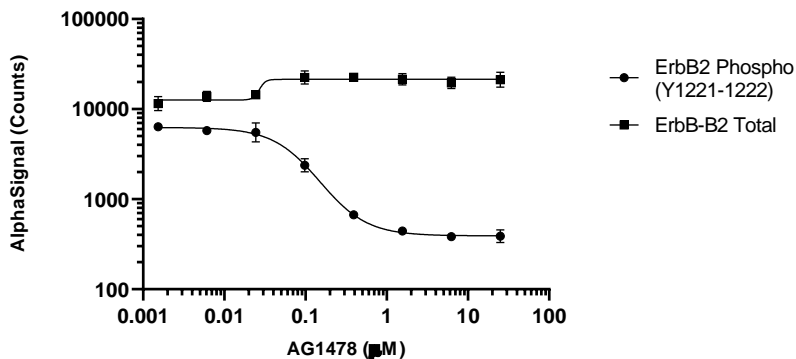
This assay kit contains antibodies which recognize invariant epitopes on Receptor tyrosine-protein kinase erbB-2. The protein detected by this kit corresponds to UniProt ID P04626. Erb B2 is also known as c-ErbB2, MLN19, Neu, HER2 and p185erbB2. These antibodies recognize ErbB2 of human, rat and mouse origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from A431 cells, cultured in a T175 flask in 10% FBS containing medium, starved overnight, then treated with 1 µg/mL EGF for 5 minutes and lysed in 4 mL of 1X *SureFire Ultra* Lysis Buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. A431 cells were plated at 40K cells per well in a 96 well plate and incubated overnight. Cells were treated with the indicated concentrations of AG1478 for 2 hours prior to being activated by 1 µg/mL EGF for 5 minutes. Cells were lysed with 1x Lysis Buffer and assayed separately for Total or Phospho ErbB2 (Y1221/1222) using respective *SureFire Ultra* kits. Equivalent to approximately 4000 cells/datapoint.



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>