

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ ErbB2 Total Assay Kit

**Products: ALSU-TEB2-A500, ALSU-TEB2-A10K, ALSU-TEB2-A50K,
ALSU-TEB2-A-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize invariant epitopes on erythroblastic leukemia viral oncogene homolog 2 (ErbB2). The protein detected by this kit corresponds to GenBank Accession NP_001005862. ErbB2 is also known as NEU, NGL, HER2, TKR1, CD340, HER-2, MLN 19 and HER-2/neu.

These antibodies recognize ErbB2 of human, rat and mouse origin. Other species should be tested on a case-by-case basis.

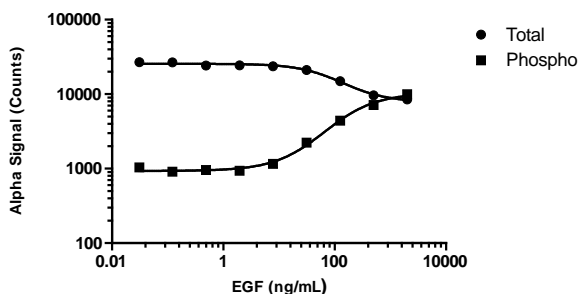
Control Lysate Information

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

Representative data

Data obtained with a 2-plate, 2-incubation protocol. Confluent A431 cells were treated with EGF at the indicated concentrations for 5 minutes, then lysed with 1X Lysis buffer. Levels of ErbB2 (Tyr1221/1222) phosphorylation and ErbB2 Total levels were assayed with respective *SureFire Ultra* assay kits.

EGF Dose Response in A431 cells - ErbB2 Ultra assays



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>