

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ p-ErbB2 (Tyr1221/1222) Assay Kit

**Products: ALSU-PEB2-A500, ALSU-PEB2-A10K, ALSU-PEB2-A50K,
ALSU-PEB2-A-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Tyr1221/1222 epitope, and a distal epitope, on v-erb-b2 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.

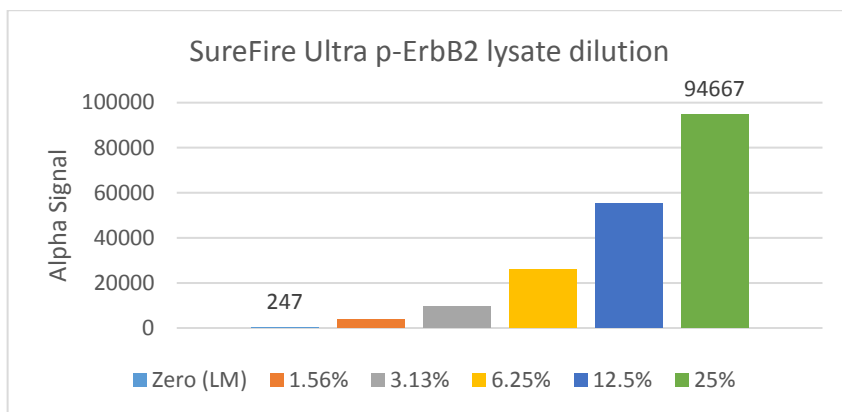
These antibodies recognize ErbB2 of human and mouse origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from A431 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, 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.



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>