

# Technical Data Sheet

Caution: Research use only.

## Alpha SureFire® Ultra™ Multiplex p-ErbB2 ( Tyr1221/1222) + Total ErbB2 Assay Kit Products: MPSU-PTEB2-K500, MPSU-PTEB2-K10K, MPSU-PTEB2-K50K, MPSU-PTEB2-K-HV

### Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Tyr1221/1222 epitope, and a distal epitope, on ErbB2, plus antibodies to measure the total levels of ErbB2. The protein detected by this kit corresponds to Entre Gene ID 2064. These antibodies recognize ErbB2 of human, rat and mouse origin. Other species should be tested on a case-by-case basis.

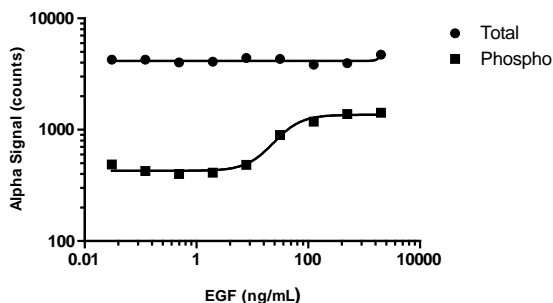
### Control Lysate Information

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 SKBR3 cells were serum starved for 2 hours, then treated with EGF at the indicated concentrations for 5min. Cells were then lysed with 1X Lysis buffer and assayed for both phospho and total ErbB2 as detailed in the MPSU manual.

#### EGF Dose Response in SKBR3 cells - ErbB2 MPSU



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>