

# Technical Data Sheet

Caution: Research use only.

## Alpha SureFire® Ultra™ Multiplex

### Terbium SureFire Ultra p-EGF Receptor (Tyr1068) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

**Products: TBSU-PEGFR-A500, TBSU-PEGFR-A10K, TBSU-PEGFR-A50K, TBSU-PEGFR-A-HV**

#### **Kit-Specificity Information**

This assay kit contains antibodies which recognize the phospho-Tyr1068 epitope, and a distal epitope, on epidermal growth factor receptor (EGFR). The protein detected by this kit corresponds to GenBank Accession NP\_005219. EGFR is also known as ERBB, HER1, mENA, ERBB1 and PIG61. These antibodies recognize EGFR of human and mouse origin. Other species should be tested on a case-by-case basis.

The antibodies in this kit are identical to those in the corresponding ALSU-PEGFR-A kits, other than the CaptSure2 tag.

#### **Control Lysate Information**

Positive Control Lysate: T175 Flasks of confluent A431 cells were serum-starved for 24 hours, then stimulated with 10µg/mL EGF for 10 minutes and lysed with 4 mL of 1X Lysis buffer.

#### **Representative data**

Control lysate samples measured in the terbium channel using an Envision plate reader.

