

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex

Terbium SureFire Ultra p-IGF Receptor β (Tyr1135/1136) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

**Products: TBSU-PIGFR-B500, TBSU-PIGFR-B10K, TBSU-PIGFR-B50K,
TBSU-PIGFR-B-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Tyr1135/1136 epitope, and a distal epitope, on the IGF1 Receptor β subunit. The protein detected by this kit correspond to GenBank Accession NP_000866. IGF1R is also known as IGFR, CD221, IGFIR, JTK13, MGC18216, MGC142170 and MGC142172. These antibodies recognize IGF1R of human, mouse and rat 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-PIGFR-B kits, other than the CaptSure2 tag.

Control Lysate Information

Positive Control Lysate: A431 cells were cultured to confluence T175 flasks. The cells were then serum-starved overnight, and treated with insulin (200 μ g/mL) for 5 minutes. The cells were lysed with 5mL of freshly prepared 1X Lysis buffer, with shaking for 10 minutes.

Representative data

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

