

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex p-IGF-1 Receptor β (Tyr1135/1136) + Total IGF-1 Receptor β Assay Kit

**Products: MPSU-PTIGFR-K500, MPSU- PTIGFR-K10K, MPSU- PTIGFR-K50K,
MPSU- PTIGFR-K-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Tyr1135/1136 epitope, and a distal epitope, on IGF-1 Receptor β for phospho IGF-1 Receptor β analysis, plus antibodies to measure the total levels of IGF-1 Receptor β .

These antibodies recognize IGF-1 Receptor β of human, mouse and rat origin. Other species should be tested on a case-by-case basis.

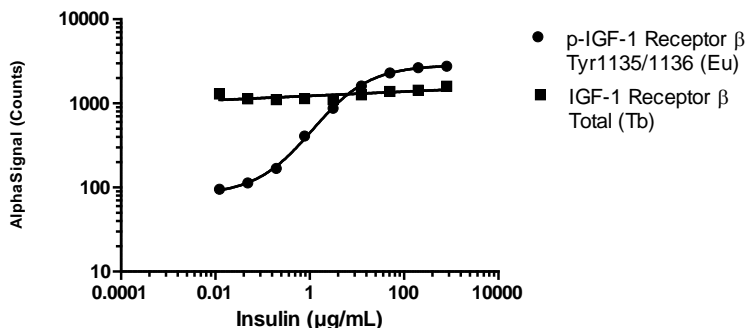
Control Lysate Information

Positive Control Lysate: Prepared from A431 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 200 μ g/mL Insulin for 5min and lysed in 4mL of 1X *SureFire Ultra* Lysis buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. A431 cells were plated at 20K cells per well in a 96 well plate, left overnight and then treated for 5 minutes with insulin in serum-free medium at the indicated concentrations. Cells were lysed with 50 μ L of 1X Lysis buffer and assayed for both phospho and total IGF-I Receptor β .

Insulin Dose Response- A431 (5min Treatment)



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>