

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex

Terbium SureFire Ultra p-Insulin Receptor β (Tyr1150/1151) Assay Kit For multiplexing with AlphaLISA SureFire Ultra Assay Kits

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

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Tyr1150/1151 epitope, and a distal epitope, on insulin receptor β subunit (INSR). The protein detected by this kit corresponds to GenBank Accession NP_000199. INSR is also known as HHF5 and CD220. These antibodies recognize INSR 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-PINR-A kits, other than the CaptSure2 tag.

Control Lysate Information

Positive Control Lysate: Confluent T175 flasks of CHO cells transfected with the insulin receptor (CHO-INSR cells) were stimulated with 2 $\mu\text{g}/\text{mL}$ insulin for 15 minutes and then lysed with 8mL of freshly prepared 1X Lysis buffer, with shaking for 10 minutes. Once reconstituted, lysates should be stored frozen in single use aliquots.

Representative data

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

