

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex p-Insulin Receptor β (Tyr1150/1151) + Total Insulin Receptor β Assay Kit

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

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Tyr1150/1151 epitope, and a distal epitope, on Insulin Receptor β for phospho Insulin Receptor β analysis, plus antibodies to measure the total levels of Insulin Receptor β .

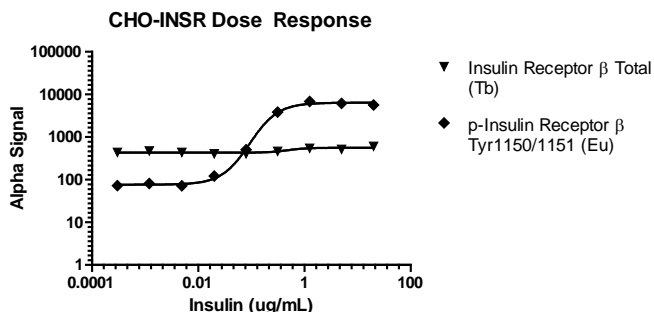
These antibodies recognize Insulin Receptor β of human, mouse and rat origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from CHO cells transfected with the insulin receptor (CHO-INSR cells), cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 2 μ g/mL Insulin and 0.1% BSA for 15 minutes and lysed in 4mL of 1X *SureFire Ultra* Lysis buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. CHO-INSR cells were plated overnight at 40K cells per well in a 96 well plate, and treated for 15 minutes with insulin in serum-free medium at the indicated concentrations. Cells were lysed with 100 μ L of 1X Lysis buffer and samples were assayed simultaneously for phospho and total Insulin Receptor β .



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>