

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ p-Insulin Receptor β (Tyr1150/1151) Assay Kit

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

Kit-Specificity Information

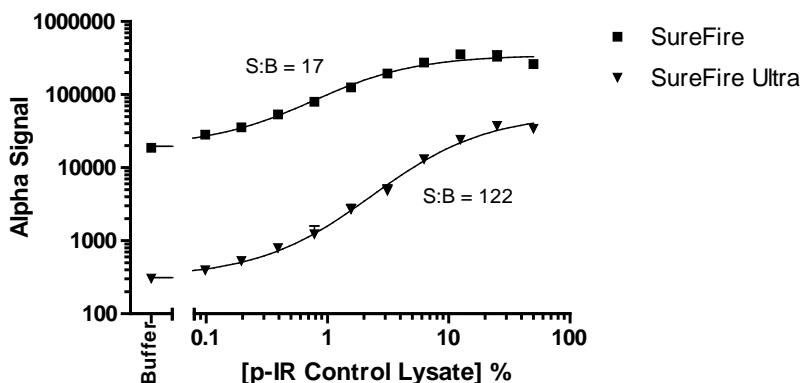
This assay kit contains antibodies which recognize the phospho-Tyr1150/1151 epitope, and a distal epitope, on insulin receptor β subunit (INSR). The protein detected by this kit corresponds to GenBank Accession NP_000199. INSR is also known as HHF5 and CD220. These antibodies recognize INSR of human, mouse and rat origin. Other species should be tested on a case-by-case basis.

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. Data shows a comparison of AlphaScreen *SureFire* and AlphaLISA *SureFire Ultra* assay performance. Control lysate was diluted in the Lysis buffer specific for the assay type.



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>