

# Technical Data Sheet

Caution: Research use only.

## AlphaLISA® SureFire® Ultra™ p-IGF-1 Receptor $\beta$ (Tyr1135/1136) Assay Kit

**Products: ALSU-PIGFR-B500, ALSU-PIGFR-B10K, ALSU-PIGFR-B50K,  
ALSU-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  $\beta$  subunit. The protein detected by this kit correspond to GenBank Accession NP\_000866. IGF1R is also known as IGFR, CD221, IGFR, 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.

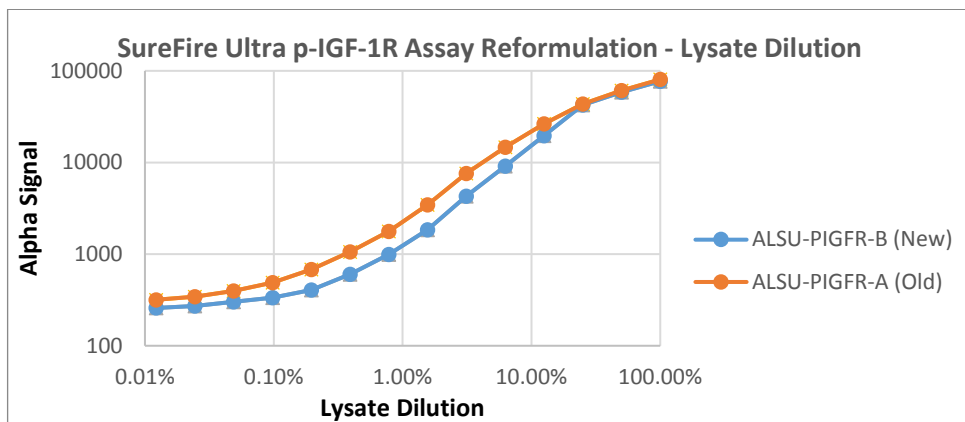
### **Control Lysate Information**

Positive Control Lysate: Prepared from A431 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 200 $\mu$ 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.

Data is shown for this new formulation and the previous old kit formulation.



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>