

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ FGF Receptor 1 Total Assay Kit

ALSU-TFGFR1-A500, ALSU-TFGFR1-A10K, ALSU-TFGFR1-A50K, ALSU-TFGFR1-A-HV

Kit-Specificity

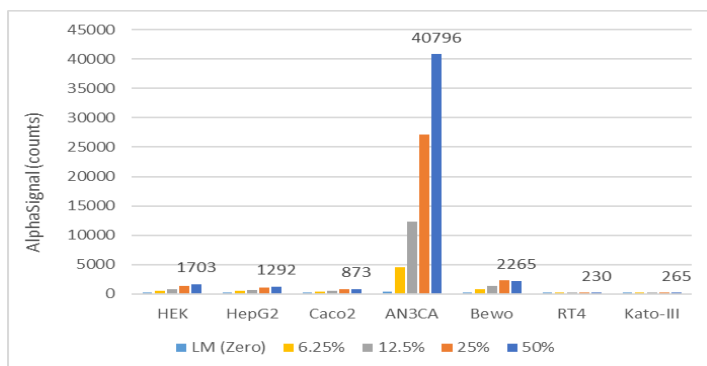
This assay kit contains antibodies which recognize invariant epitopes on Fibroblast Growth Receptor 1 (FGFR1). The protein detected by this kit corresponds to UniProt ID P11362. FGFR1 is also known as bFGF-R-1, FLT-2, N-sam and c-Fgr. These antibodies recognize FGF Receptor 1 of human origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from AN3-CA cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 0.5 mM Pervanadate for 10 minutes, and lysed in 5 mL 1X *SureFire Ultra* Lysis Buffer.

Representative Data

Data obtained with a 2 plate, 2 incubation protocol. Data obtained for Total FGFR-1 from screening various lysates prepared from cells lysed in *SureFire Ultra* Lysis Buffer. The '50%' point corresponds to approximately 5,000 cells/datapoint for Caco2, Bewo and Kato-III cells up to 25,000 cells/datapoint for AN3CA cells. HEK, HepG2 and RT4 cells/datapoint were approximately 10,000 at the '50%' level.



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>.