

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ FGF Receptor 3 Total Assay Kit

ALSU-TGFR3-A500, ALSU-TGFR3-A10K, ALSU-TGFR3-A50K, ALSU-TGFR3-A-HV

Kit-Specificity

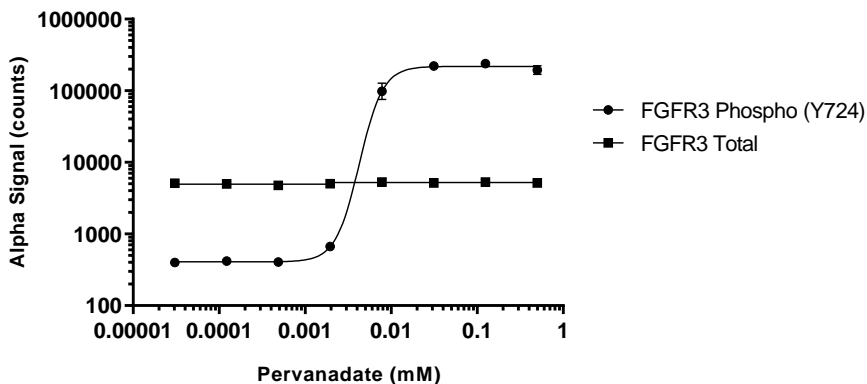
This assay kit contains antibodies which recognize two invariant epitopes on FGF Receptor 3. The protein detected by this kit corresponds to UniProt ID P22607. These antibodies recognize FGF Receptor 3 of human origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from HepG2 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 10 mL 1X *SureFire Ultra* Lysis Buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. HepG2 cells were plated at 40K cells per well in a 96 well plate and incubated overnight. Cells were then treated with the Pervanadate at the indicated concentrations for 10 minutes. Cells were lysed with 1x Lysis Buffer and assayed separately for Total and Phospho (Tyr724) FGF Receptor 3, using respective *SureFire Ultra* kits. Equivalent to approximately 4000 cells/datapoint.



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>.