

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ p-FGF Receptor 2 (Tyr653/654) Assay Kit

**Products: ALSU-PFGFR2-A500, ALSU-PFGFR2-A10K, ALSU-PFGFR2-A50K,
ALSU-PFGFR2-A-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Tyr653/654 epitope, and a distal epitope on FGF Receptor 2. The protein detected by this kit corresponds to Entrez-Gene ID 2263. These antibodies recognize FGF Receptor 2 of human origin. Other species should be tested on a case-by-case basis.

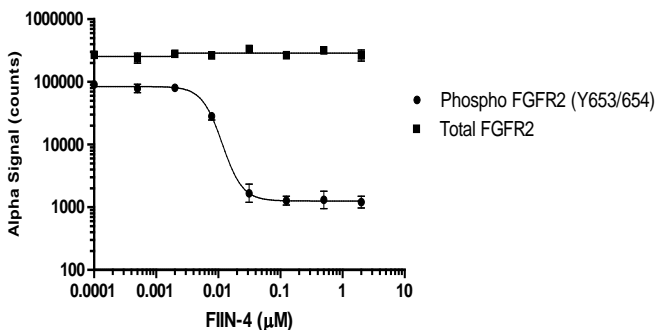
Control Lysate Information

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. KATO-III cells were plated overnight at 40K cell/well, then inhibited with the indicated concentrations of FIIN-4 for 16 hours. Cells were lysed with 100 μ L of 1X Lysis Buffer and assayed separately for total and phospho FGF Receptor 2, using respective kits.

FIIN-4 Treatment Of KATO-III Cells: pFGFR2 (Y653/654)



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>.