

Technical Data Sheet

Caution: Research use only.

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

**Products: ALSU-TFGFR3-A500, ALSU-TFGFR3-A10K, ALSU-TFGFR3-A50K,
ALSU-TFGFR3-A-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize total levels of FGF Receptor 3. The protein detected by this kit corresponds to Entrez-Gene ID 2261. These antibodies recognize FGF Receptor 3 of human origin. Other species should be tested on a case-by-case basis.

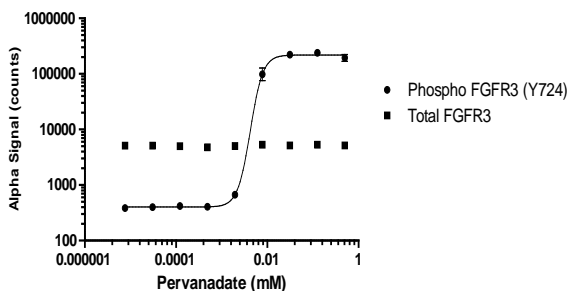
Control Lysate Information

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 overnight at 40K cell/well, then inhibited with the indicated concentrations of Pervanadate for 10 minutes. Cells were lysed with 120 µL of 1X Lysis Buffer and assayed separately for total and phospho FGF Receptor 3, using respective kits.

Pervanadate Dose Response In HepG2 Cells: pFGFR3 (Y724)



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>.