

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ p-FGF Receptor 3 (Tyr724) Assay Kit

**Products: ALSU-PFGFR3-A500, ALSU-PFGFR3-A10K, ALSU-PFGFR3-A50K,
ALSU-PFGFR3-A-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Tyr724 epitope, and a distal epitope on FGF Receptor 3. The protein detected by this kit corresponds to Entrez-Gene ID 2261. These antibodies recognize FGF Receptor 3 of human and mouse origin. Other species should be tested on a case-by-case basis.

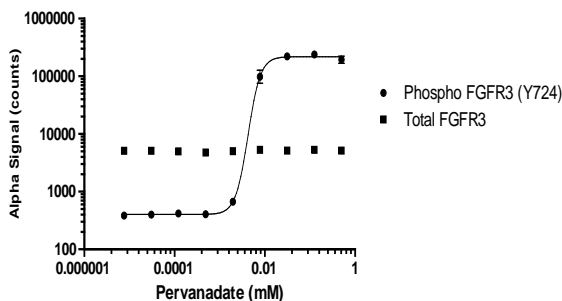
Control Lysate Information

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 μ L 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>.