

AlphaLISA® SureFire® Ultra™

Human FGF Receptor 1 Total Detection Kit

Product number: ALSU-TFGFR1-A500, ALSU-TFGFR1-A10K,
 ALSU-TFGFR1-A50K, ALSU-TFGFR1-A-HV



Kit specificity:

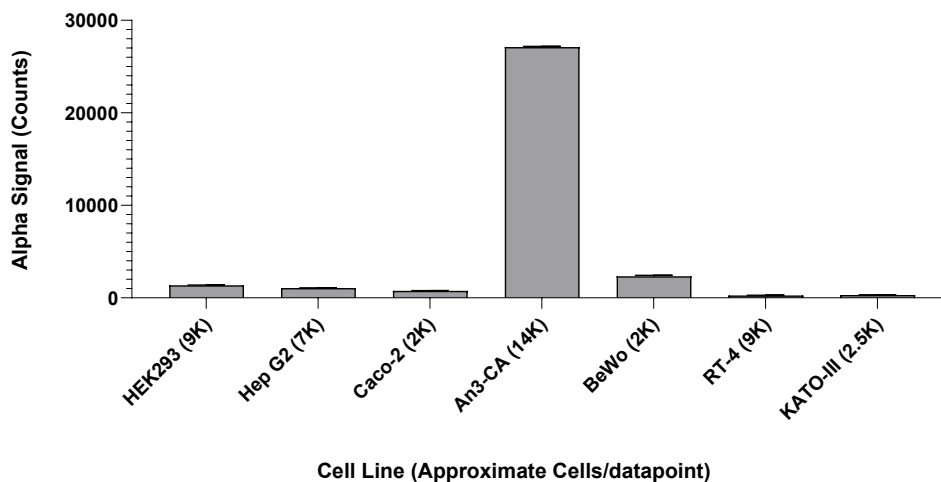
This assay kit contains antibodies which recognize distinct epitopes on FGFR1. The protein detected by this kit corresponds to UniProt ID P11362. FGFR1 is also known as bFGF-R-1 and FLT-2. These antibodies recognize FGF Receptor 1 of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

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

Representative data

Data obtained from measurement of FGF Receptor 1 Total in various cell types lysed with Lysis Buffer. Approximate number of cells/datapoint is indicated for the various cell lines.



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF"