

AlphaLISA® SureFire® Ultra™

Human p-VEGF Receptor 2 (Tyr1214) Detection Kit

Product number: ALSU-PVGFR-C500, ALSU-PVGFR-C10K,
ALSU-PVGFR-C50K, ALSU-PVGFR-C-HV



Kit specificity:

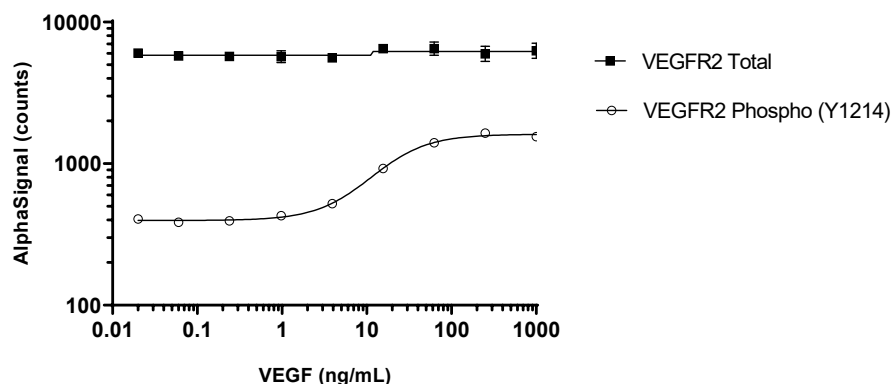
This assay kit contains antibodies which recognize the phospho-Tyr1214 epitope and a distal epitope, on VEGFR2. The protein detected by this kit corresponds to UniProt ID P35968. VEGFR2 is also known as Vascular endothelial growth factor receptor 2, FLK1 and KDR. These antibodies recognize VEGFR2 of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

Positive Control Lysate: Prepared from HUVEC cells, cultured to confluence in T175 flasks in 2% FBS containing medium. Cells were serum starved for 2 hours in 0.5% FBS containing media, treated with 100 ng/mL rh-VEGF for 5 minutes and lysed with 4 mL of Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. HUVEC cells were plated at 20K cells/well in a 96 well plate and incubated overnight. Cells were starved overnight prior to treatment with rh-VEGF at the indicated concentrations for 5 minutes. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Tyr1214) and Total VEGFR2 using respective *SureFire Ultra* kits. Equivalent to approximately 2,000 cells/datapoint.



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"