

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ VEGF Receptor 2 Total Assay Kit

ALSU-TVGFR-A500, ALSU-TVGFR-A10K, ALSU-TVGFR-A50K, ALSU-PVGFR-A-HV

Kit-Specificity

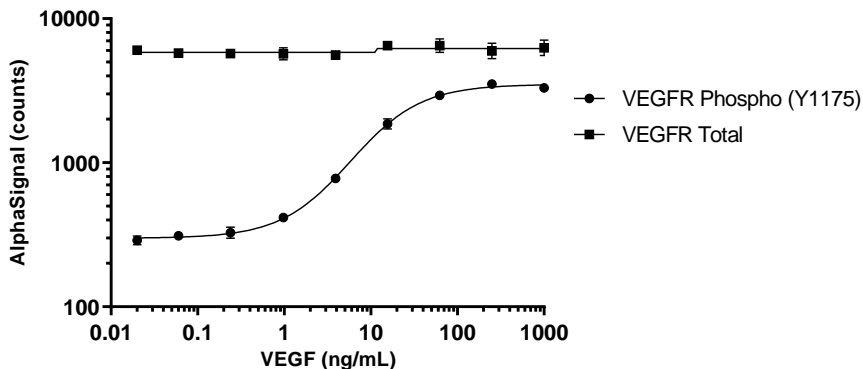
This assay kit contains antibodies which recognize the two invariant epitopes on vascular endothelial growth factor receptor 2 (VEGFR2). The protein detected by this kit corresponds to UniProt ID P35968. VEGFR2 is also known as FLK1, CD309 and KDR. These antibodies recognize VEGFR2 of human origin. Other 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 medium, starved for 2 hours, then treated with 100 ng/mL rh-VEGF for 5 minutes and lysed in 4 mL of 1X *Surefire Ultra* Lysis Buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. HUVEC cells were plated at 20K cell per 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 1x Lysis Buffer and assayed separately for Total or Phospho (Tyr1175) VEGFR2 using respective *SureFire Ultra* kits. Equivalent to approximately 2000 cells/datapoint.



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>