

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ p-VEGF Receptor 2 (Tyr1175) Assay Kit

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

Kit-Specificity Information

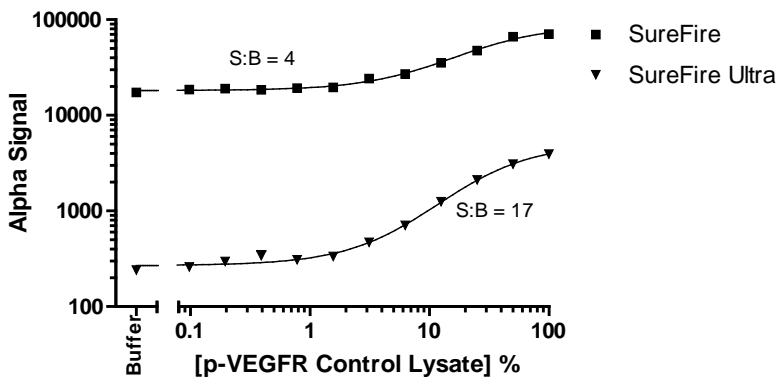
This assay kit contains antibodies which recognize the phospho-Tyr1175 epitope, and a distal epitope, on vascular endothelial growth factor receptor 2 (VEGFR2). The protein detected by this kit corresponds to GenBank Accession NP_002244. VEGFR2 is also known as FLK1, CD309, VEGFR, and KDR. These antibodies 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 100ng/mL rhVEGF for 5 minutes and lysed in 4mL of 1X *Surefire Ultra* lysis buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. Data shows a comparison of AlphaScreen *SureFire* and AlphaLISA *SureFire Ultra* assay performance. Control lysate was diluted in the Lysis buffer specific for the assay type.



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>