

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ p-VEGF Receptor 2 (Tyr1214) Assay Kit

**Products: ALSU-PVGFR-C500, ALSU-PVGFR-C10K, ALSU-PVGFR-C50K,
ALSU-PVGFR-C-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Tyr1214 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.

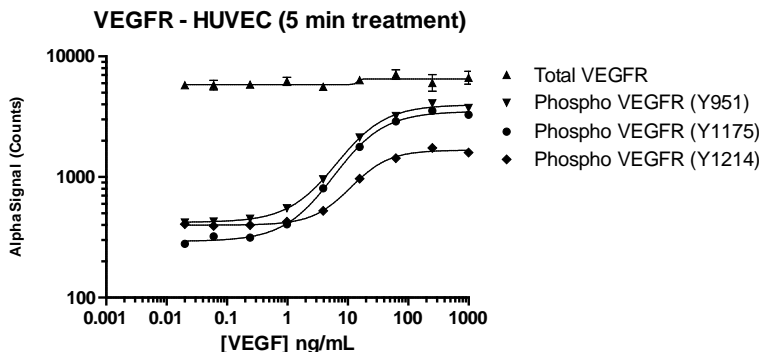
The human Tyr1214 site is equivalent to the mouse Tyr1212 site.

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. HUVEC cells were grown to confluence, starved for 3 hours, and stimulated with rh-VEGF for 5 min at 37C.



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>