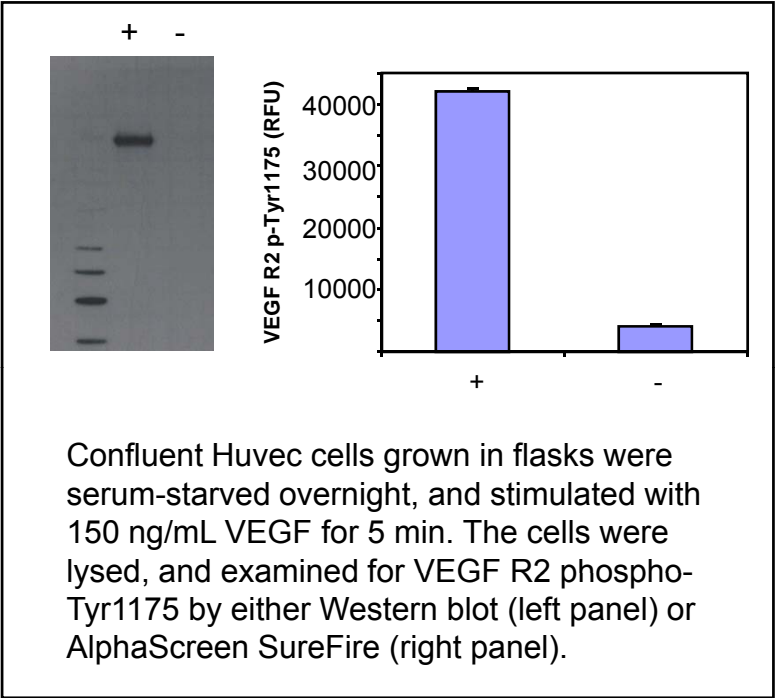
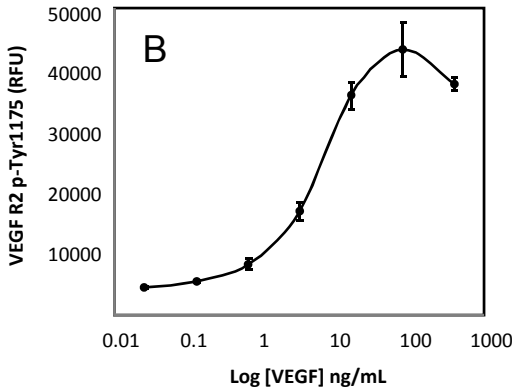
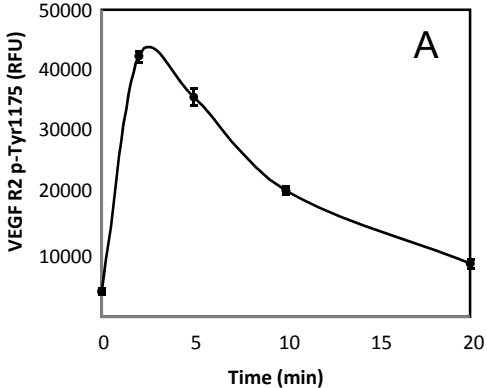


TGRVRS500TK
TGRVRS500
TGRVRS10K
TGRVRS50K

# AlphaScreen® SureFire® phospho-VEGF receptor 2 (Tyr1175) Kit

Vascular endothelial growth factor receptor 2 (VEGF R2) is a major receptor for VEGF-mediated signaling in endothelial cells. Upon ligand binding, VEGF R2 undergoes autophosphorylation in the kinase insert domain (Tyr951/996) and in the tyrosine kinase catalytic domain (Tyr1054/1059) to become activated. Activation of the receptor leads to rapid recruitment of adaptor proteins. The phosphorylation of Tyr1175 mediates binding with the p85 subunit of PI-3 kinase, PLCγ, and Shb.



HUVEC cells were seeded at 15,000 cells per well into 96-well microplates, and grown to confluence over 3 days. The cells starved overnight in serum-free media. The following day the media was removed, and the cells were treated with (A) 1000 ng/mL VEGF for various lengths of time, or (B) treated with various concentrations of VEGF for 10 min. The agonist was removed, and the cells were lysed with 50 μL per well of freshly prepared 1X Lysis buffer, with gentle shaking for 10 min. The lysate was subsequently analysed for VEGF R2 phospho-Tyr1175 using the standard AlphaScreen SureFire protocol.