

# Technical Data Sheet

Caution: Research use only.

## Alpha SureFire® Ultra™ Multiplex

### Terbium SureFire Ultra p-VEGF Receptor 2 (Tyr1175) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

**Products: TBSU-PVGFR-A500, TBSU-PVGFR-A10K, TBSU-PVGFR-A50K, TBSU-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.

The antibodies in this kit are identical to those in the corresponding ALSU-PVGFR-A kits, other than the CaptSure2 tag.

#### **Control Lysate Information**

Positive Control Lysate: HUVEC cells were grown to confluence in T175 flasks, and then serum-starved overnight. The following day the media was removed, and the cells were treated with 100ng/mL VEGF for 10 minutes. The agonist was removed, and the cells were lysed with 4mL of freshly prepared 1X Lysis buffer, with shaking for 10 minutes.

#### **Representative data**

Control lysate samples measured in the terbium channel using an Envision plate reader.

