

AlphaLISA® SureFire® Ultra™

Human p-TIE2 (Tyr992) Detection Kit

Product number: ALSU-PTIE2-A500, ALSU-PTIE2-A10K
 ALSU-PTIE2-A50K, ALSU-PTIE2-A-HV



Kit specificity:

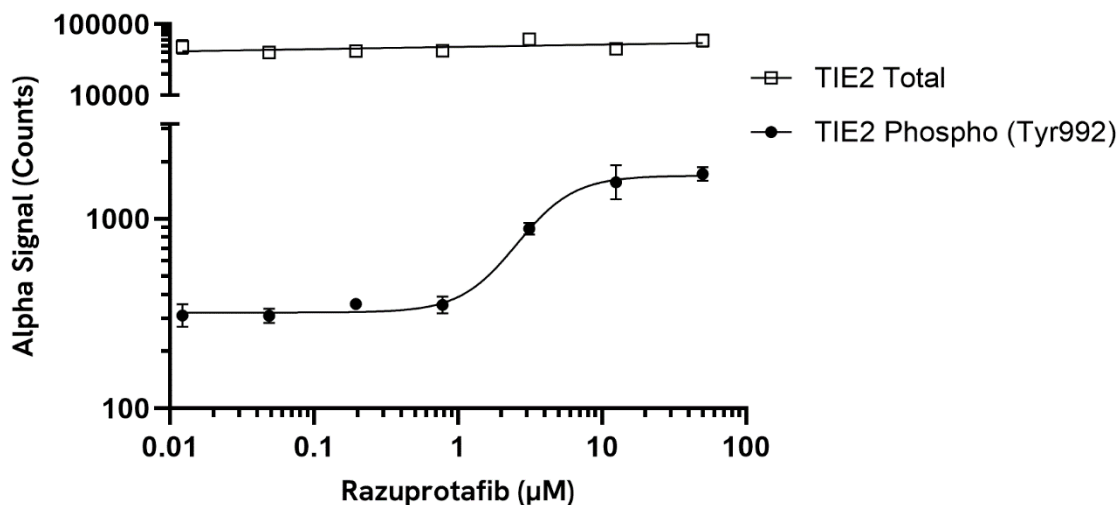
This assay kit contains antibodies which recognize the phospho-Tyr992 epitope and a distal epitope on TIE2. The protein detected by this kit corresponds to UniProt ID Q06807. TIE2 is also known as Angiopoietin-1 receptor and TEK. These antibodies recognize TIE2 of human origin. Other species should be tested on a case-by case basis.

Control lysate information:

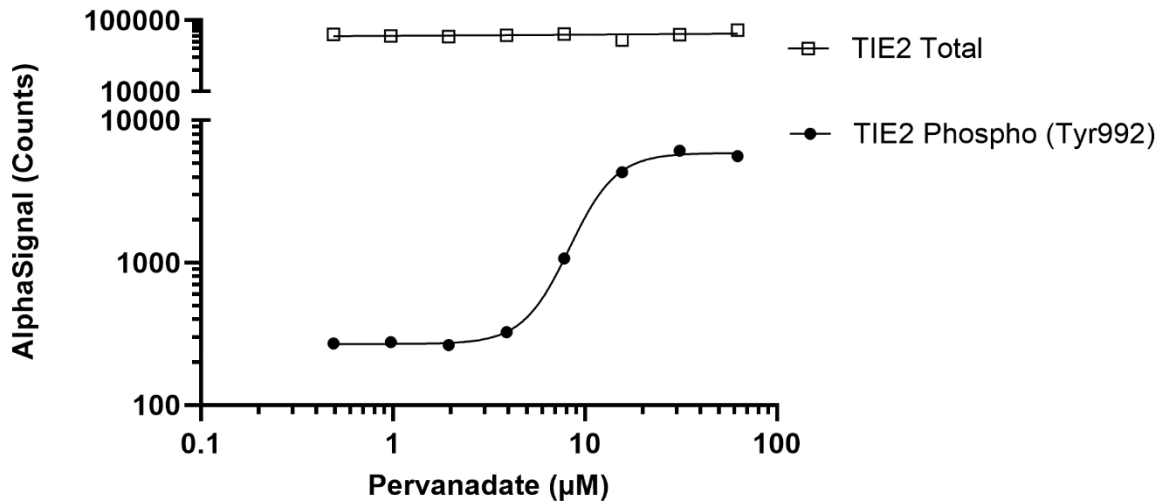
Positive Control Lysate: Prepared from HUVEC cells cultured to confluence in a T175 cm flask in FBS containing medium, then treated with 250 µM Pervanadate for 30 minutes and lysed with 4 mL of Lysis Buffer.

Representative data:

Data obtained with a 2-plate, 2-incubation protocol. HUVEC cells were seeded at 20K cells/well in a 96 well plate and incubated overnight. Cells were starved for 2 hours and then treated with Razuprotafib at the indicated concentrations for 15 minutes. Cells were lysed and assayed separately for Phospho (Tyr992) and Total TIE2 using respective SureFire Ultra kits. Equivalent to approximately 4,000 cells/datapoint.



Data obtained with a 2-plate, 2-incubation protocol. HUVEC cells were seeded at 20K cells/well in a 96 well plate and incubated for 48 hours. Cells were treated with Pervanadate at the indicated concentrations for 30 minutes. Cells were lysed and assayed separately for Phospho (Tyr992) and Total TIE2 using respective *SureFire Ultra* kits. Equivalent to approximately 8,000 cells/datapoint.



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF"

