Research use only. Not for use in diagnostic procedures.

AlphaLISA® SureFire® Ultra™ Human FLT3 Total Detection Kit

Product number:

ALSU-TFLT3-A500, ALSU-TFLT3-A10K,

ALSU-TFLT3-A50K, ALSU-TFLT3-A-HV



Kit specificity:

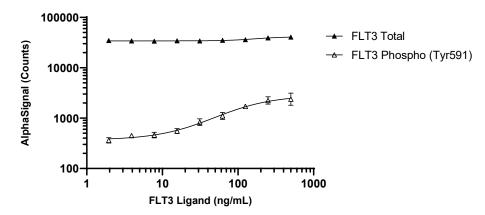
This assay kit contains antibodies which recognize distinct epitopes on FLT3. The protein detected by this kit corresponds to UniProt ID P36888. FLT3 is also known as FL cytokine receptor, FLK-2 and STK-1. These antibodies recognize FLT3 of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

Positive Control Lysate: Prepared from THP 1 cells cultured in 10% FBS containing media. Cells were harvested, washed with HBSS + 0.1% BSA and adjusted to $10 \times 10^{\circ}$ cells/mL. Cells were treated with 500 ng/mL FLT3L for 15 minutes and lysed with 5X Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. THP-1 cells were washed, resuspended in HBSS + 0.1% BSA and seeded at 250K cells/well in a 96 well plate. Cells were treated with FLT3L at the indicated concentrations for 10 minutes. Cells were lysed with the addition of 5X Lysis Buffer and assayed separately for Phospho (Tyr591) and Total FLT3 using respective *SureFire Ultra* kits. Equivalent to approximately 20,000 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"





Revvity, Inc. 940 Winter Street Waltham, MA 02451 USA (800) 762-4000 www.revvity.com For a complete listing of our global offices, visit <u>www.revvity.com</u> Copyright ©2023, Revvity, Inc. All rights reserved.