

Research use only. Not for use in diagnostic procedures.

AlphaLISA® SureFire® Ultra™

Human p-TFEB (Ser122) Detection Kit

Product number:

ALSU-PTFEB-A500, ALSU-PTFEB-A10K,

ALSU-PTFEB-A50K, ALSU-PTFEB-A-HV



Kit specificity:

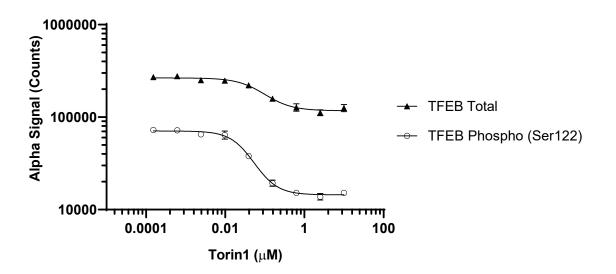
This assay kit contains antibodies which recognize the phospho-Ser122 epitope and a distal epitope on TFEB. The protein detected by this kit corresponds to UniProt ID P19484. TFEB is also known as TCFEB and transcription factor EB. These antibodies recognize TFEB of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

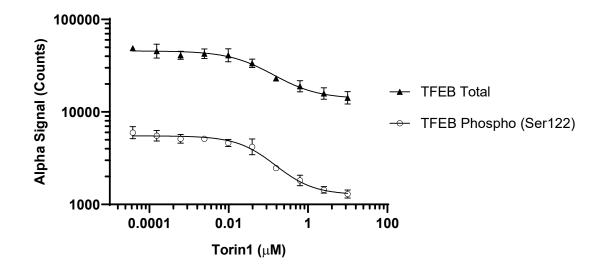
Positive Control Lysate: Prepared from A431 cells, cultured to confluency in T175 flasks in 10% FBS containing medium and lysed with 10 mL of Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. A431 cells were plated at 60K cells/well in a 96 well plate and incubated overnight. Cells were treated with Torin1 at the indicated concentrations for 3 hours. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Ser122) and Total TFEB using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint.



Data obtained with a 2-plate, 2-incubation protocol. HepG2 cells were plated at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with Torin1 at the indicated concentrations for 3 hours. Cells were lysed with Lysis Buffer and assayed separately for TFEB Phospho (Ser122) and TFEB Total using respective *SureFire Ultra* kits. Equivalent to approximately 2,600 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.