

AlphaLISA® SureFire® Ultra™

Human and Mouse GAPDH Total Detection Kit

Product number: ALSU-TGAPD-B500, ALSU-TGAPD-B10K,
ALSU-TGAPD-B50K, ALSU-TGAPD-B-HV



Kit specificity:

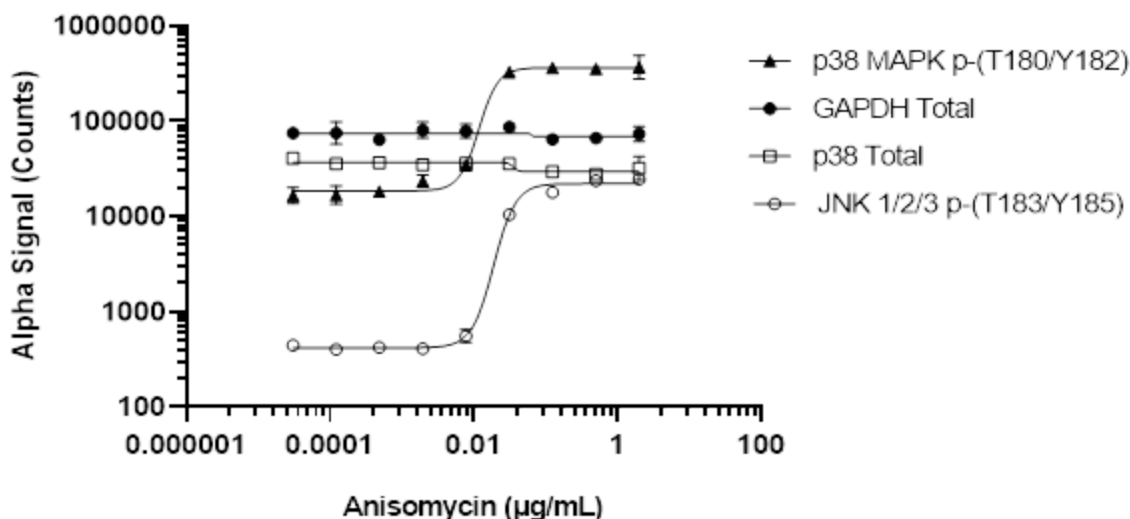
This assay kit contains antibodies which recognize distinct epitopes on GAPDH. The protein detected by this kit corresponds to UniProt ID P04406. GAPDH is also known as Peptidyl-cysteine S-nitrosylase GAPDH. These antibodies recognize GAPDH of human and mouse origin. Other species should be tested on a case-by-case basis.

Control lysate information:

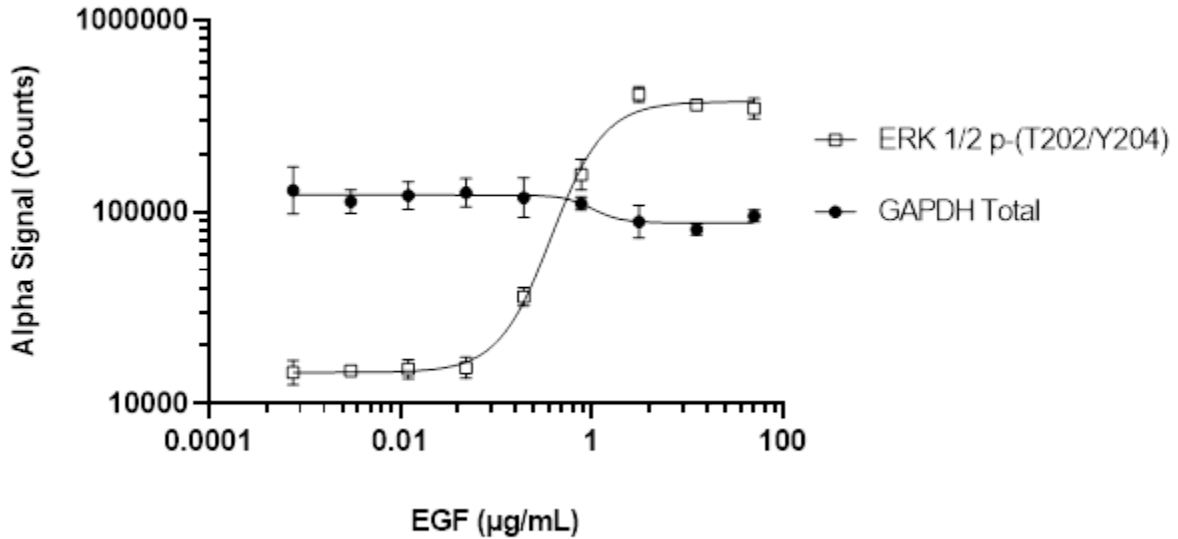
Positive Control Lysate: Prepared from HeLa cells, cultured to confluence in T175 flasks in 10% FBS containing medium, lysed with 40 mL of Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. NIH3T3 cells were seeded at 30K cells/well in a 96 well plate and incubated overnight. Cells were serum deprived for 2 hours then treated with Anisomycin at the indicated concentrations for 30 minutes. Cells were lysed with Lysis Buffer and assayed separately for Phospho-(Thr180/Tyr182) p38 MAPK, Total p38, Phospho-(Thr183/Tyr185) JNK 1/2/3 and Total GAPDH using respective *SureFire Ultra* kits. Equivalent to approximately 3,000 cells/datapoint for Phospho and Total p38, Phospho JNK 1/2/3 and 1,000 cells/datapoint for Total GAPDH.



Data obtained with a 2-plate, 2-incubation protocol. HeLa cells were seeded at 50K cells/well in a 96 well plate and incubated overnight. Cells were serum deprived for 2 hours and then treated with EGF at the indicated concentrations for 10 minutes. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Thr202/Tyr204) ERK and Total GAPDH using respective *SureFire Ultra* kits. Equivalent to approximately 2,500 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"

