

AlphaLISA® SureFire® Ultra™ Terbium

Human and Mouse Cofilin Total Detection Kit

Product number: TBSU-TCOF-A500, TBSU-TCOF-A10K
TBSU-TCOF-A50K, TBSU-TCOF-A-HV



Kit specificity:

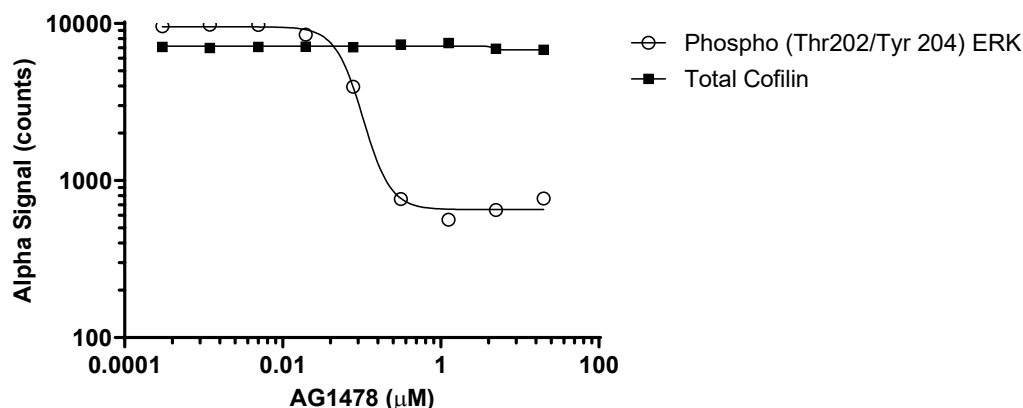
This assay kit contains antibodies which recognize distinct epitopes on Cofilin. The protein detected by this kit corresponds to UniProt ID P23528. Cofilin is also known as CFL. These antibodies recognize Cofilin of human and mouse origin. Other species should be tested on a case-by-case basis.

Control lysate information:

Positive Control Lysate: Prepared from HeLa cells, cultured to confluence in T175 flasks in 10% FBS containing media and lysed with 12 mL of Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. HeLa cells were seeded at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with AG1478 at the indicated concentrations for 2 hours and then treated with 1 ng/mL of EGF for 10 minutes. Cells were lysed with Lysis buffer and assayed for Phospho (Thr202/Tyr204) ERK and Total Cofilin in the same well of an assay plate using the SureFire Ultra (Phospho (Thr202/Tyr204) ERK kit paired with the Terbium SureFire Ultra Total Cofilin kit.



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"