

## Alpha SureFire<sup>®</sup> Ultra<sup>™</sup> Multiplex

# Human and Mouse p-Cofilin (Ser3) + Total Cofilin Detection Kit

**Product number:** 

MPSU-PTCOF-K500, MPSU-PTCOF-K10K,

MPSU-PTCOF-K50K, MPSU-PTCOF-K-HV



#### Kit specificity:

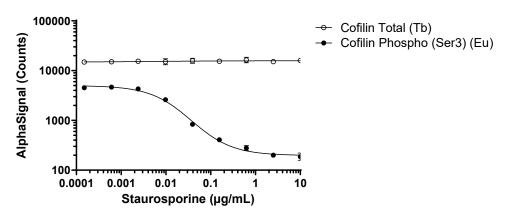
This assay kit contains antibodies which recognize the phospho-Ser3 epitope and a distal epitope on Cofilin, plus antibodies to measure the total levels of Cofilin. The protein detected by this kit corresponds to UniProt ID P23528. Cofilin is also known as CFL1. These antibodies recognize cofilin of 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 20K cells per well in a 96 well plate and grown overnight. Cells were treated with Staurosporine at the indicated concentrations for 60 minutes. Cells were lysed with Lysis buffer and assayed for Phospho (Ser3) and Total Cofilin in the same wells of an assay plate using the *SureFire Ultra* Multiplex kit. Equivalent to 1,000 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.

VersionTGRSU288.1