

# Technical Data Sheet

Caution: Research use only.

## Alpha SureFire® Ultra™ Multiplex p-Cofilin (Ser3) + Total Cofilin Assay Kit

**Products: MPSU-PTCOF-K500, MPSU-PTCOF-K10K, MPSU-PTCOF-K50K,  
MPSU-PTCOF-K-HV**

### Kit-Specificity Information

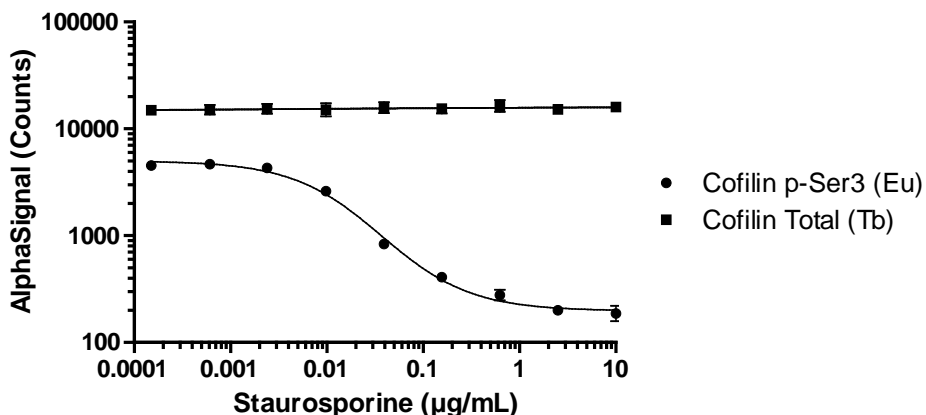
This assay kit contains antibodies which recognize the phospho-Ser3 epitope, and a distal epitope, on cofilin for phospho cofilin analysis, plus antibodies to measure the total levels of cofilin. The protein detected by this kit corresponds to GenBank Accession NP\_005498.1. Also known as CFL, CFL1. These antibodies recognize cofilin of human, mouse and rat 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 and lysed in 12ml of 1X *SureFire Ultra* Lysis Buffer.

### Representative data

Data obtained with a 2-plate, 2-incubation protocol. HeLa cells were plated overnight at 20K cells per well in a 96 well plate, and then treated with staurosporine at the indicated concentrations for 60 minutes. Cells were lysed with 200µL of 1X Lysis buffer and samples were assayed for both total and phosphor (Ser3) Cofilin, as detailed in the MPSU manual.



For more information and detailed protocols, please refer to the generic AlphaLISA *SureFire Ultra* Assay Kits Manual, available from <http://www.perkinelmer.com/category/alpha-surefire-kits>