

# Technical Data Sheet

Caution: Research use only.

## AlphaLISA® SureFire® Ultra™ Cofilin Total Assay Kit

**Products: ALSU-TCOF-A500, ALSU-TCOF-A10K, ALSU-TCOF-A50K,  
ALSU-TCOF-A-HV**

### **Kit-Specificity Information**

This assay kit contains antibodies which recognize two invariant epitopes on cofilin. The protein detected by this kit corresponds to GenBank Accession NP\_005498.1. Also known as CFL, CFL1.

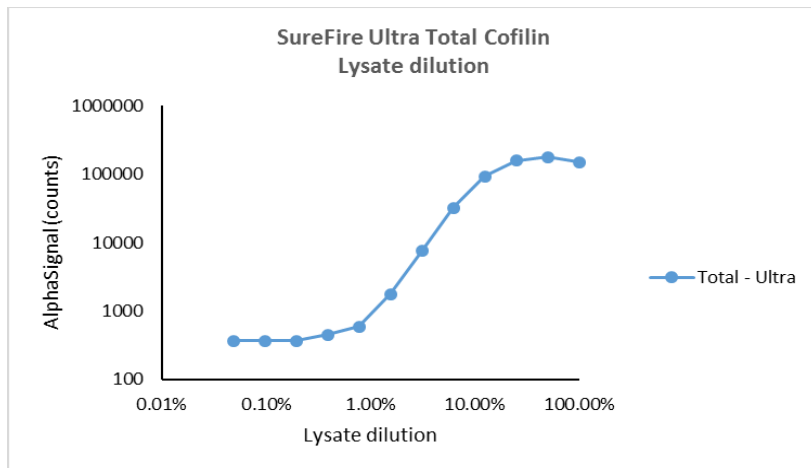
These antibodies recognize cofilin of human, mouse and rat origin. Other species should be tested on a case-by-case basis.

### **Control Lysate Information**

Positive Control Lysate: Prepared from HeLa cells, cultured to confluence in T175 flasks 10% FBS containing medium and lysed in 12ml 1X *SureFire Ultra* Lysis Buffer.

### **Representative data**

Data obtained with a 2-plate, 2-incubation protocol.



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>