

# Technical Data Sheet

Caution: Research use only.

## Alpha SureFire® Ultra™ Multiplex p-eIF2α (Ser51) + Total eIF2α Assay Kit

**Products: MPSU-PTEIF2-K500, MPSU-PTEIF2-K10K, MPSU-PTEIF2-K50K,  
MPSU-PTEIF2-K-HV**

### Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Ser51 epitope, and a distal epitope, on eukaryotic translation initiation factor 2 subunit 1 (eIF2α), plus antibodies to measure the total levels of eIF2α. The protein detected by this kit corresponds to GenBank Accession NP\_004085. eIF2α is also known as EIF2A, EIF-2A, EIF-2alpha and EIF2S1.

These antibodies recognize eIF2α of human and mouse origin. Other species should be tested on a case-by-case basis.

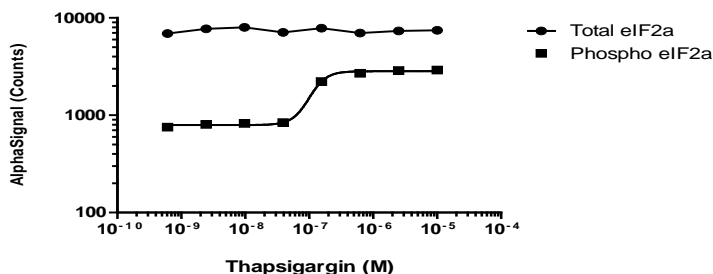
### Control Lysate Information

Positive Control Lysate: Prepared from C2C12 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 1μM Thapsigargin for 30min and lysed in 4mL of 1X *SureFire Ultra* Lysis buffer.

### Representative data

Data obtained with a 2-plate, 2-incubation protocol. Confluent C2C12 cells were treated with Thapsigargin at the indicated concentrations for 60min, then lysed with 1X Lysis buffer. Samples were assayed for both Phospho and Total eIF2α as detailed in the MPSU manual.

#### **Thapsigargin Dose Response in C2C12 cells - eIF2a MPSU**



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>