

# Technical Data Sheet

## AlphaLISA® SureFire® Ultra™ p-eIF2α (Ser51) Assay Kit

ALSU-PEIF2-B500, ALSU-PEIF2-B10K, ALSU-PEIF2-B50K, ALSU-PEIF2-B-HV

### Kit-Specificity

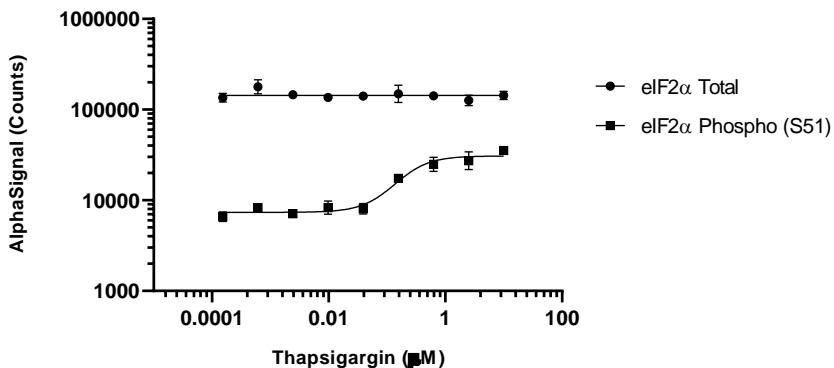
This assay kit contains antibodies which recognize the phospho-Ser51 epitope, and a distal epitope, on eukaryotic translation initiation factor 2 subunit 1 (eIF2α). The protein detected by this kit corresponds to UniProt ID Q9BY44. These antibodies recognize eIF2α of human and mouse origin. Other species 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 100 nM Thapsigargin for 30 minutes and lysed in 4 mL of 1X *SureFire Ultra* Lysis Buffer.

### Representative Data

Data obtained with a 2-plate, 2-incubation protocol. C2C12 cells were plated at 20K cells per well in a 96 well plate and incubated overnight. Cells were then treated with Thapsigargin at the indicated concentrations for 60 minutes. Cells were lysed with 1x Lysis Buffer and assayed separately for Total or Phospho (Ser51) eIF2α using respective *SureFire Ultra* kits. Equivalent to approximately 2000 cells/datapoint.



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>