

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex

Terbium SureFire Ultra p-eIF2 α (Ser51) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

**Products: TBSU-PEIF2-B500, TBSU-PEIF2-B10K, TBSU-PEIF2-B50K,
TBSU-PEIF2-B-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 (eIF2 α). The protein detected by this kit corresponds to GenBank Accession NP_004085. eIF2 α is also known as EIF2A, EIF-2A, EIF-2alpha and EIF2S1. These antibodies recognize eIF2 α of human, rat and mouse origin. Other species should be tested on a case-by-case basis.

The antibodies in this kit are identical to those in the corresponding ALSU-PEIF2-B kits, other than the CaptSure2 tag.

Control Lysate Information

Positive Control Lysate: Prepared from C2C12 cells, cultured to confluence in T175 flasks in 10% FBS containing medium for 3 days, then treated with thapsigargin (1 μ M) for 30min and lysed in 4mL of 1X SureFire Ultra Lysis buffer.

Representative data

Control lysate samples measured in the terbium channel using an Envision plate reader.

