

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex

Terbium SureFire Ultra p-eIF4E (Ser209) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

**Products: TBSU-PEIF4-A500, TBSU-PEIF4-A10K, TBSU-PEIF4-A50K,
TBSU-PEIF4-A-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Ser209 epitope, and a distal epitope, on eukaryotic translation initiation factor 4E (eIF4E). The protein detected by this kit corresponds to GenBank Accession NP_001124150. eIF4E is also known as CBP, EIF4F, EIF4E1, EIF4EL1 and MGC111573. These antibodies recognize eIF4E of human, mouse and rat 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-PEIF4-A kits, other than the CaptSure2 tag.

Control Lysate Information

Positive Control Lysate: Prepared from HEK293 cells, cultured to confluence in T175 flasks in 10% FBS containing medium for 3 days, then treated with EGF (200ng/mL) for 30min and lysed in 12mL of 1X SureFire Ultra Lysis buffer.

Representative data

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

