

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex

Terbium SureFire Ultra p-CREB (Ser133) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

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

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Ser133 epitope, and a distal epitope, on cAMP responsive element binding protein 1 (CREB). The protein detected by this kit corresponds to GenBank Accession NP_004370. CREB is also known as MGC9284 or CREB1. These antibodies recognize CREB 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-PCREB-A kits, other than the CaptSure2 tag.

Control Lysate Information

Positive Control Lysate: Prepared from A431 cells, cultured to confluence in T175 flasks in 10% FBS containing medium for 3 days, then treated with EGF (500ng/mL) for 10min and lysed in 8mL of 1X SureFire Ultra Lysis buffer.

Representative data

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

