

# Technical Data Sheet

Caution: Research use only.

## AlphaLISA® SureFire® Ultra™ CREB Total Assay Kit

**Products: ALSU-TCREB-A500, ALSU-TCREB-A10K, ALSU-TCREB-A50K,  
ALSU-TCREB-A-HV**

### **Kit-Specificity Information**

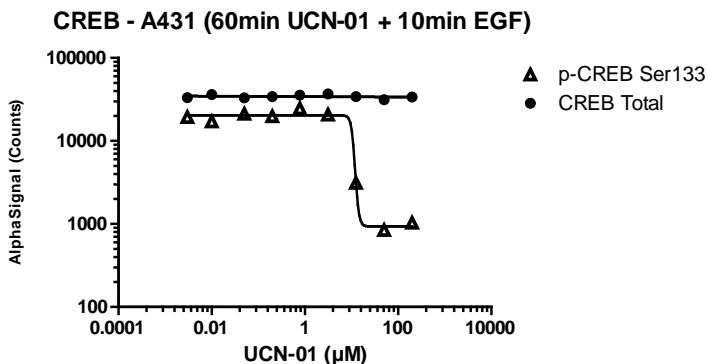
This assay kit contains antibodies which recognize two invariant epitopes on cAMP responsive element binding protein 1 (CREB). The protein detected by this kit corresponds to GenBank Accession NP\_004370. CREB is also known as MGC9284 or CREB1. These antibodies recognize CREB of human, mouse and rat origin. Other species should be tested on a case-by-case basis.

### **Control Lysate Information**

Positive Control Lysate: Prepared from A431 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, starved for 60min, then treated with 500ng/mL EGF for 10min and lysed in 8mL of 1X *SureFire Ultra* Lysis buffer.

### **Representative data**

Data obtained with a 2-plate, 2-incubation protocol. A431 cells were treated with the protein kinase inhibitor UCN-01 at the indicated concentrations for 60min, then with EGF (200ng/mL) for 10min. Cells were then lysed with 1X Lysis Buffer. Samples were separately assayed for p-CREB (Ser133) and CREB Total, using respective *SureFire Ultra* kits.



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>