

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ p-CREB (Ser133) Assay Kit

ALSU-PCREB-A500, ALSU-PCREB-A10K, ALSU-PCREB-A50K, ALSU-PCREB-A-HV

Kit-Specificity

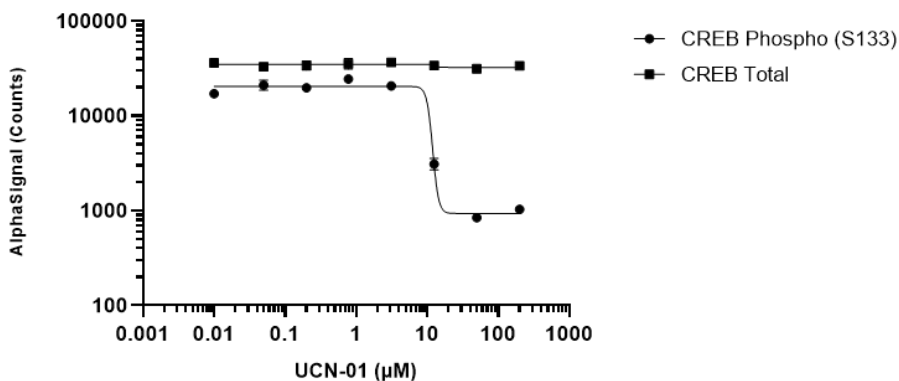
This assay kit contains antibodies which recognize the phospho Ser-133 epitope, and a distal epitope on CREB. The protein detected by the kit corresponds to UniProt ID P16220. CREB is also known as cAMP-responsive element-binding protein 1. These antibodies recognize CREB of human, mouse and rat origin. 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 60 minutes, then treated with 500 ng/mL EGF for 10 minutes and lysed in 8 mL 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 60 minutes at 37°C, then with 200 ng/mL EGF for 10 minutes. Cells were then lysed with 1X Lysis Buffer. Samples were separately assayed for Total or Phospho (Ser133) CREB, with respective *SureFire Ultra* kits. Equivalent to approximately 12,000 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>