

# Technical Data Sheet

Caution: Research use only.

## Alpha SureFire® Ultra™ p-CHK1 (Ser345) Assay Kit

**Products: ALSU-PCHK1-A500, ALSU-PCHK1-A10K, ALSU-PCHK1-A50K,  
ALSU-PCHK1-A-HV**

### Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Ser345 epitope, and a distal epitope, on CHK1. The protein detected by this kit corresponds to GenBank Accession NP\_001107594. CHK1 is also known as CHEK1. These antibodies recognize CHK1 of human and mouse origin. Other species should be tested on a case-by-case basis.

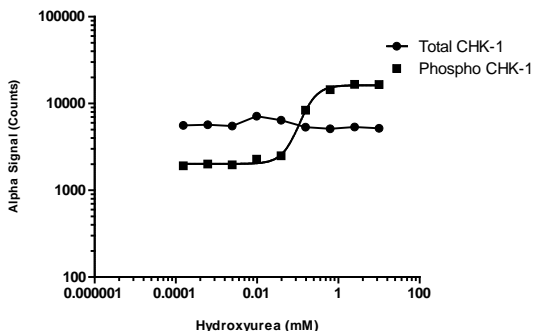
### Control Lysate Information

Positive Control Lysate: Prepared from HEK293 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 5mM hydroxyurea for 2 hours, and lysed in 4mL of 1X *SureFire Ultra* Lysis buffer.

### Representative data

Data obtained with a 2-plate, 2-incubation protocol. HEK293 cells were plated overnight at 40K cells per well in a 96 well plate, and treated for 2 hours in serum-containing medium with hydroxyurea at the indicated concentrations. Cells were lysed with 100µL of 1X Lysis buffer and assayed separately for phospho and total CHK1 with respective *SureFire Ultra* assay kits.

Hydroxyurea Dose Response - HEK 293 cells



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>