

# Technical Data Sheet

## AlphaLISA® SureFire® Ultra™ Histone H3 Total Assay Kit

**ALSU-THISH3-A500, ALSU-THISH3-A10K, ALSU-THISH3-A50K, ALSU-THISH3-A-HV**

### **Kit Specificity**

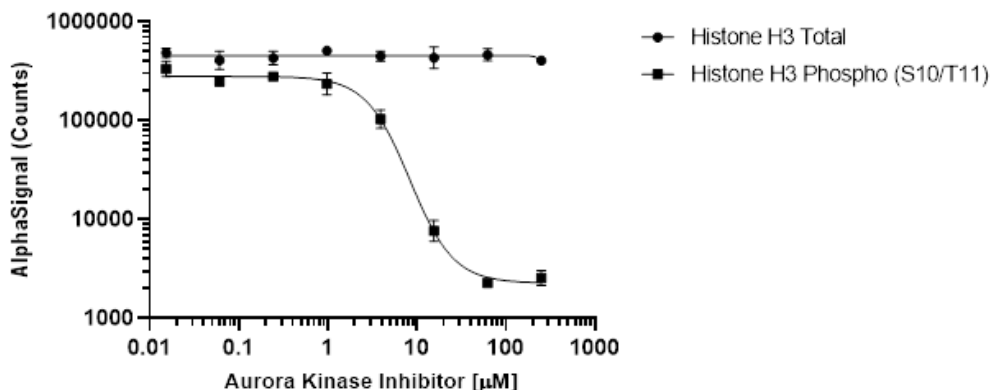
This assay kit contains antibodies which recognize invariant epitopes on Histone H3. The protein detected by this kit corresponds to UniProt ID P68431. These antibodies recognize Histone H3 of human and mouse origin. Other species should be tested on a case-by-case basis.

### **Control Lysate Information**

Positive Control Lysate: Prepared from U2OS cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 10  $\mu$ M nocodazole for 16 hours and lysed with the equivalent of 2000 mL of 1X *SureFire Ultra* Lysis Buffer C.

### **Representative Data**

Data obtained with a 2-plate, 2-incubation protocol. U2OS cells were plated at 5K cells per well in a 96 well plate and incubated overnight. Cells were then treated with Aurora Kinase Inhibitor at the indicated concentrations for 30 minutes before the addition of 100 nM nocodazole for a further 16 hours. Cells were lysed with 1X *SureFire Ultra* Lysis Buffer C and assayed separately for Total and Phospho (Ser10/Thr11) Histone H3, using respective *SureFire Ultra* assay kits. Equivalent to approximately 100 cells/datapoint. Nb: Histone H3 is expressed at high levels in mammalian cells so extensive titration of cells/lysates is recommended to ensure operation in the linear range of the assay.



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>