

AlphaLISA® SureFire® Ultra™

Human and Mouse Histone H3 Total Detection Kit

Product number: ALSU-THISH3-A500, ALSU-THISH3-A10K,
 ALSU-THISH3-A50K, ALSU-THISH3-A-HV



Kit specificity:

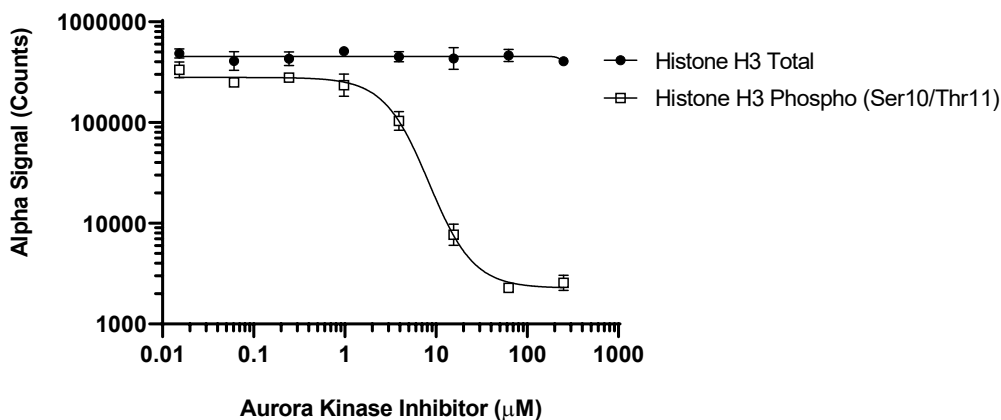
This assay kit contains antibodies which recognize distinct epitopes on Histone H3. The protein detected by this kit corresponds to UniProt ID P68431. Histone H3 is also known as Histone H3.1. These antibodies recognize Histone H3 of human and mouse origin. Other species should be tested on a case-by-case 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 μ M Nocodazole for 16 hours, lysed with 120 mL of Lysis Buffer C.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. U2OS cells were seeded at 5K cells/well in a 96 well plate and incubated overnight. Cells were treated with Aurora Kinase Inhibitor at the indicated concentrations for 30 minutes, then treated with 100 nM Nocodazole for 16 hours. Cells were lysed with Lysis Buffer C and assayed separately for Phospho (Ser10/Thr11) and Total Histone H3 using respective *SureFire Ultra* kits. Equivalent to approximately 100 cells/datapoint.



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF"