

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ Histone H2AX Total Assay Kit

ALSU-TH2X-A500, ALSU-TH2X-A10K, ALSU-TH2X-A50K, ALSU-TH2X-A-HV

Kit Specificity

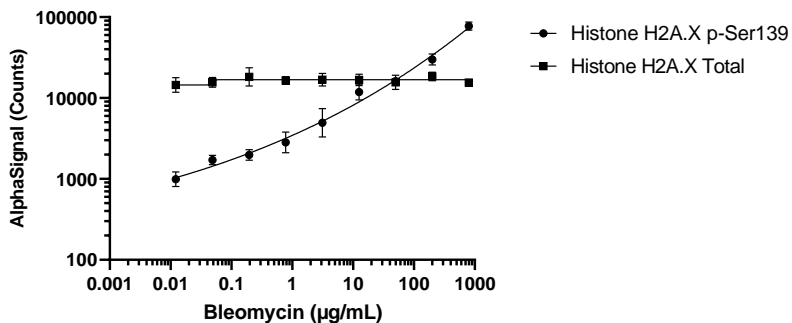
This assay kit contains antibodies which recognize invariant epitopes on Histone H2AX. The protein detected by this kit corresponds to UniProt ID P16104. Histone H2AX is also known as Histone H2A.X, H2A histone family member X, H2A.X, H2AX and H2A.X variant histone. These antibodies recognize Histone H2A.X of human origin. Other species should be tested on a case-by-case basis but given the high sequence homology, mouse reactivity is expected.

Control Lysate Information

Positive Control Lysate: Prepared from HEK293 cells cultured in T175 flasks to confluence in medium containing 10% FBS and lysed in 20mL 1X *SureFire Ultra Lysis* buffer containing 1X *SureFire Ultra Activation* buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. MCF7 cells were seeded at 10K cells/well overnight in a 96 well plate in MEM + 10% FBS. Cells were treated for 2 hrs with a dose range of bleomycin in growth media. Cells were lysed with 1X *SureFire Ultra Lysis* buffer supplemented with Activation Buffer and assayed for Phospho (Ser139) and Total Histone H2AX using respective *SureFire Ultra* kits. Equivalent to approximately 500 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>.