

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ ZAP-70 Total Assay Kit

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

Kit-Specificity

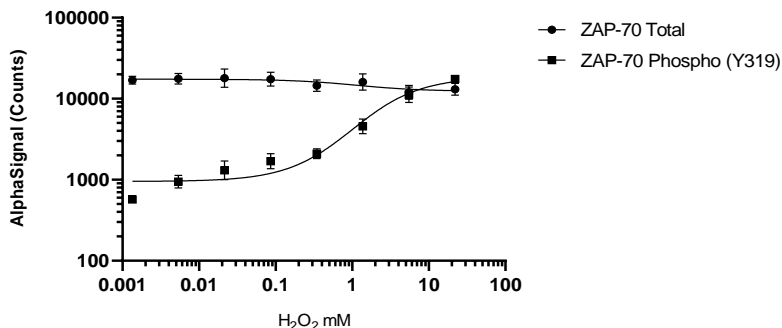
This assay kit contains antibodies that specifically recognize two invariant epitopes on ZAP-70. The protein detected by this kit corresponds to UniProt ID P43403. ZAP-70 is also known as Tyrosine-protein kinase ZAP-70 and Syk-related tyrosine kinase. These antibodies recognize ZAP-70 of human and mouse origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from Jurkat cells cultured in a T175 flask to 1×10^6 cells/mL in RPMI containing 10% FBS. Cells were treated with 11 mM H_2O_2 and 50 nM Calyculin A for 10 minutes and lysed at 3.2×10^6 cells/mL with 1X *SureFire Ultra* Lysis Buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. Jurkat cells were plated at 2×10^6 cells/mL in a 96 well plate in HBSS. Cells were then treated with H_2O_2 at the indicated concentrations for 10 minutes. Cells were then lysed with 5x Lysis Buffer at a final concentration of 0.8×10^6 cells/mL. Cell lysates were further diluted and assayed for Total or Phospho (Tyr319) Zap-70 using respective *SureFire Ultra* Kits. Equivalent to approximately 1000 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>