

Technical Data Sheet

Caution: Research use only.

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

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

Kit-Specificity Information

This assay kit contains antibodies that specifically recognize two invariant epitopes on ZAP-70. The human protein detected by this kit corresponds to Entrez Gene ID 7535. These antibodies recognize ZAP-70 of human and mouse origin. Other species should be tested on a case-by-case basis.

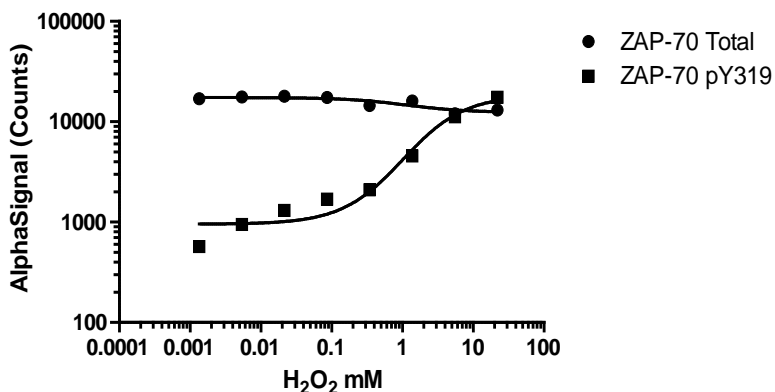
Control Lysate Information

Positive Control Lysate: Produced from Jurkat cells cultured in a T175 flask to 10^6 cells/mL in medium with 10%FBS, then treated with 11 mM H_2O_2 for 10 mins, 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 (0.2×10^6 cells/well) stimulated for 10min with H_2O_2 . Cells were lysed with 5X Lysis buffer in culture medium. Data shows signal from the equivalent of 1000 cells.

Jurkat Dose Reponse H2O2



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>