

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ KAP1 Total Assay Kit

**Products: ALSU-TKAP1-A500, ALSU-TKAP1-A10K, ALSU-TKAP1-A50K,
ALSU-TKAP1-A-HV**

Kit-Specificity Information

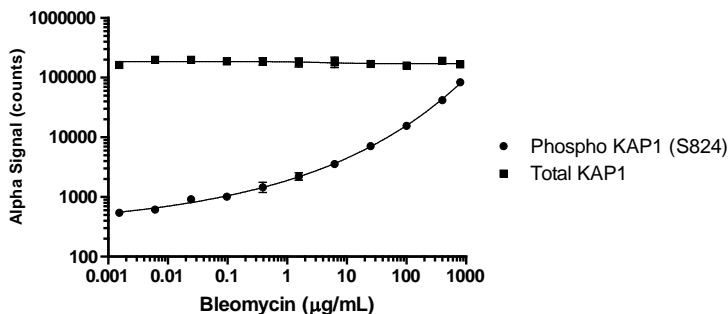
This assay kit contains antibodies which recognize invariant epitopes on KAP1. The protein detected by this kit corresponds to Entrez-Gene ID 10155. KAP1 is also known as TRIM28, TF1B and RNF26. These antibodies recognize KAP1 of human, mouse and rat origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from HEK293 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 50 µg/mL Bleomycin for 4 hours, and lysed in 8 mL 1X *SureFire Ultra* Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. HeLa cells were plated overnight at 20 K cells/well, then inhibited with the indicated concentrations of Bleomycin for 2 hours. Cells were lysed with 100 µL of 1X Lysis Buffer and assayed separately for phospho and total KAP1, using respective kits.



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>.