

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ Ras Total Assay Kit

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

Kit-Specificity

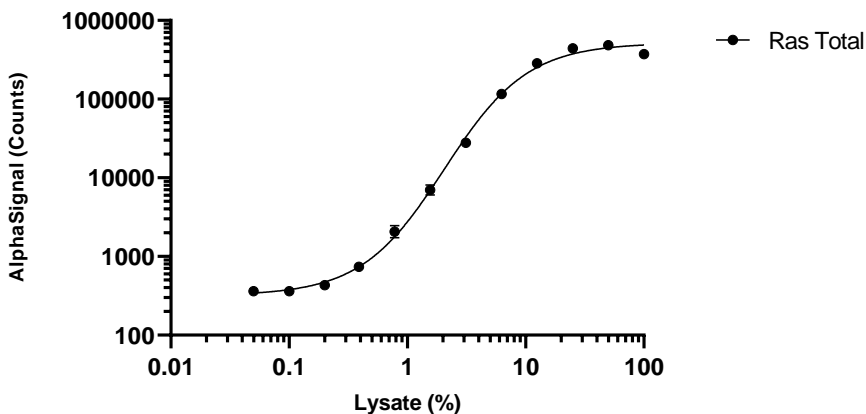
This assay kit contains antibodies which recognize invariant epitopes on Ras. The protein detected by this kit corresponds to UniProt ID P01116 (K-Ras), P01112 (H-Ras) and P01111 (N-Ras). These antibodies recognize Ras of human, rat and mouse origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from A431 cells, cultured to 80% confluence in T175 flasks in 10% FBS containing medium, and lysed in 8 mL of 1X *SureFire Ultra* Lysis Buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. A431 Positive Control Lysate was serially diluted in 1x Lysis Buffer and assayed for Total Ras using the *SureFire Ultra* kit. 100% lysate is equivalent to approximately 6000 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>