

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ GAPDH Total Assay Kit

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

Kit-Specificity

This assay kit contains antibodies which recognize invariant epitopes on glyceraldehyde-3-phosphate dehydrogenase (GAPDH). The protein detected by this kit corresponds to UniProt ID P04406. This kit recognizes GAPDH of human origin. It does not react with mouse GAPDH. Other species should be tested on a case-by-case basis.

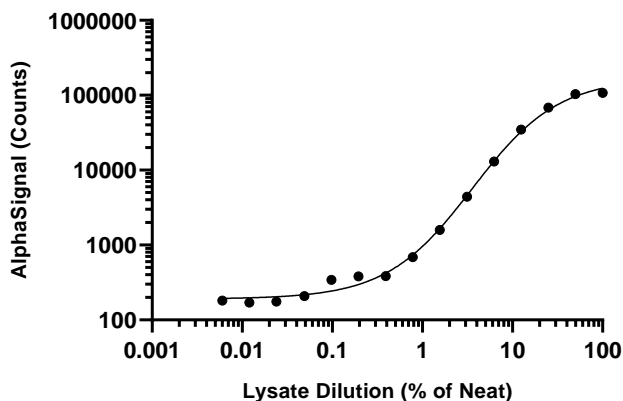
Control Lysate Information

Positive Control Lysate: Prepared from Hela cells, cultured to confluence in T75 flasks in 10% FBS containing medium, then lysed in 20mL of 1X *SureFire Ultra* Lysis Buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. Control lysate was serially diluted in 1x Lysis Buffer and assayed using the *SureFire Ultra* kit. 100% Lysate is equivalent to approximately 2000 cells/datapoint.

Note: Client should carry out a dilution of their samples to ensure they are assaying on the linear part of the curve. GAPDH is often expressed at a high concentration in mammalian cells



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>