

# Technical Data Sheet

## AlphaLISA® SureFire® Ultra™ GAPDH Total (Mouse/Human reactive) Assay Kit

ALSU-TGAPD-B500, ALSU-TGAPD-B10K, ALSU-TGAPD-B50K, ALSU-TGAPD-B-HV

### Kit-Specificity Information

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. These antibodies recognize GAPDH of human and mouse origin. Other species should be tested on a case-by-case basis.

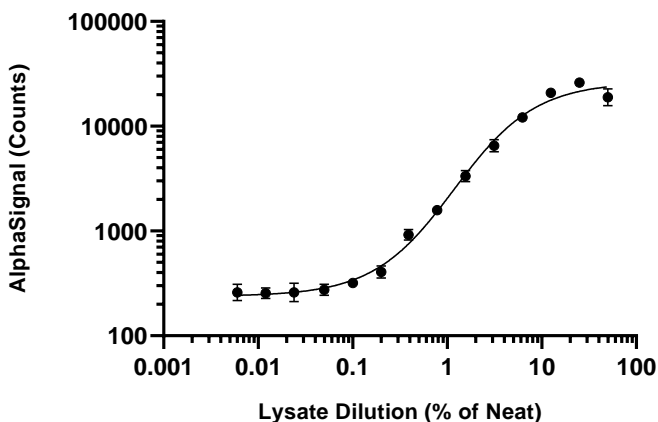
### Control Lysate Information

Positive Control Lysate: Prepared from HeLa cells, cultured to confluence in T75 flasks in 10% FBS containing medium, then lysed in 20 mL 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>