

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ p-GSK-3β (Ser9) Assay Kit

**Products: ALSU-PGS3B-A500, ALSU-PGS3B-A10K, ALSU-PGS3B-A50K,
ALSU-PGS3B-A-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Ser9 epitope, and a distal epitope, on glycogen synthase kinase 3 beta (GSK-3β). The protein detected by this kit corresponds to GenBank Accession NP_001139628. These antibodies recognize GSK-3β of human, mouse and rat origin. Other species should be tested on a case-by-case basis.

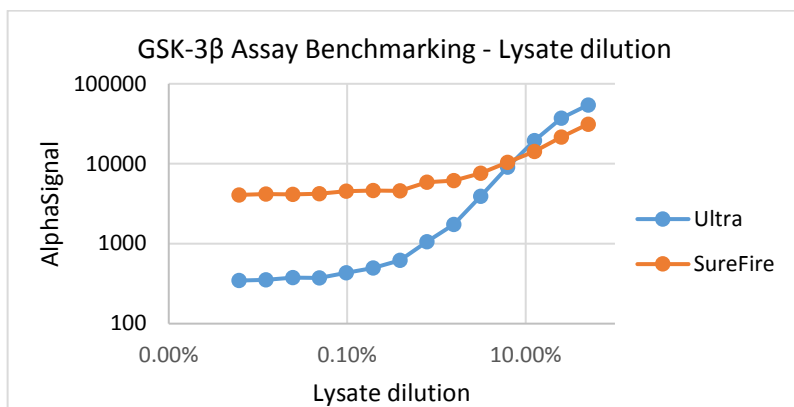
Control Lysate Information

Positive Control Lysate: Prepared from MCF7 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 500μg/mL Insulin for 30min and lysed in 20mL of 1X *SureFire Ultra* Lysis buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. Data shows a comparison of AlphaScreen *SureFire* and AlphaLISA *SureFire Ultra* assay performance. Control lysate was diluted in the Lysis buffer specific for the assay type.

S:B *SureFire Ultra* = 129; S:B *SureFire* = 8.



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>