

Technical Data Sheet

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

ALSU-PGS3B-A500, ALSU-PGS3B-A10K, ALSU-PGS3B-A50K, ALSU-PGS3B-A-HV

Kit-Specificity

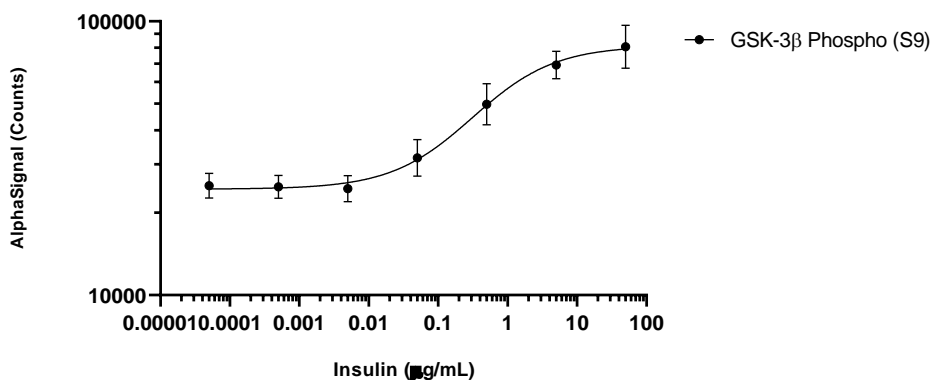
This assay kit contains antibodies which recognize the phospho-Ser9 epitope, and a distal epitope, on glycogen synthase kinase 3 beta (GSK-3β). The protein detected by this kit corresponds to UniProt ID P49841. These antibodies recognize GSK-3β of human, mouse and rat origin. Other species should be 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 500 μg/mL insulin for 30 minutes and lysed in 20 mL of 1X *SureFire Ultra* Lysis Buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. NIH/3T3 cells were plated at 40K cells per well in a 96 well plate and incubated overnight. Cells were serum starved for 2 hours and then treated with insulin at the indicated concentrations for 15 minutes. Cells were lysed with 1x Lysis Buffer and then assayed for Phospho GSK-3β (Ser9) by *SureFire Ultra*. Equivalent to approximately 4000 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>