

AlphaLISA® SureFire® Ultra™

Human and Mouse GSK3β Total Detection Kit

Product number: ALSU-TGS3B-A500, ALSU-TGS3B-A10K,
ALSU-TGS3B-A50K, ALSU-TGS3B-A-HV



Kit specificity:

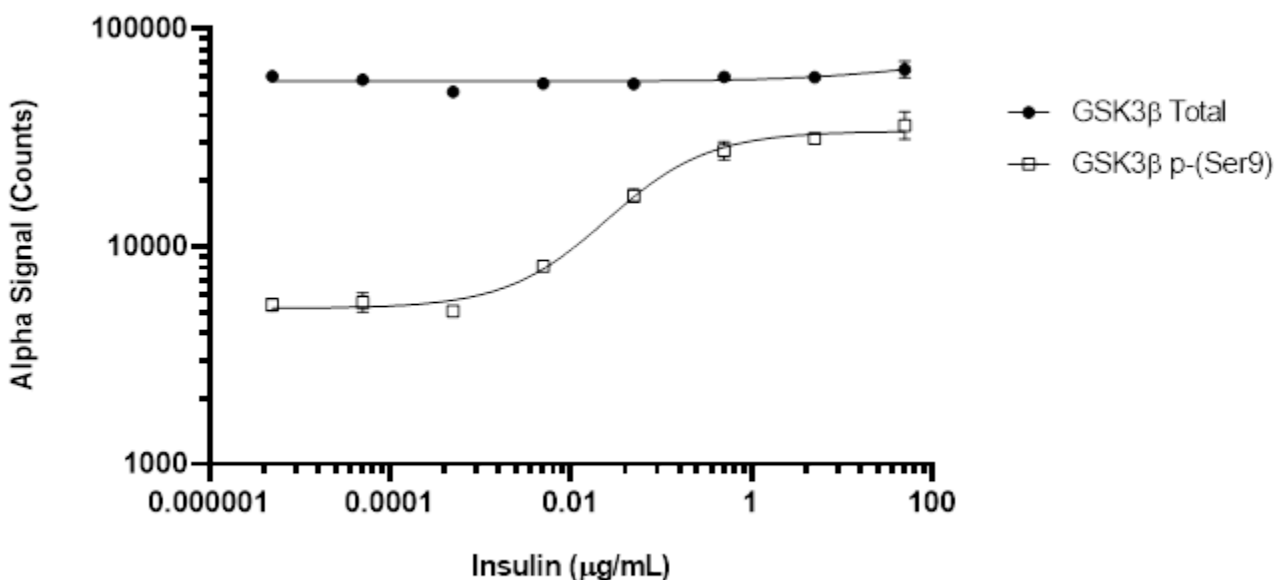
This assay kit contains antibodies which recognize distinct epitopes on GSK3β. The protein detected by this kit corresponds to UniProt ID P49841. GSK3β is also known as Serine/threonine-protein kinase GSK3β. These antibodies recognize GSK3β of human and mouse origin. Other species should be tested on a case-by-case basis.

Control lysate information:

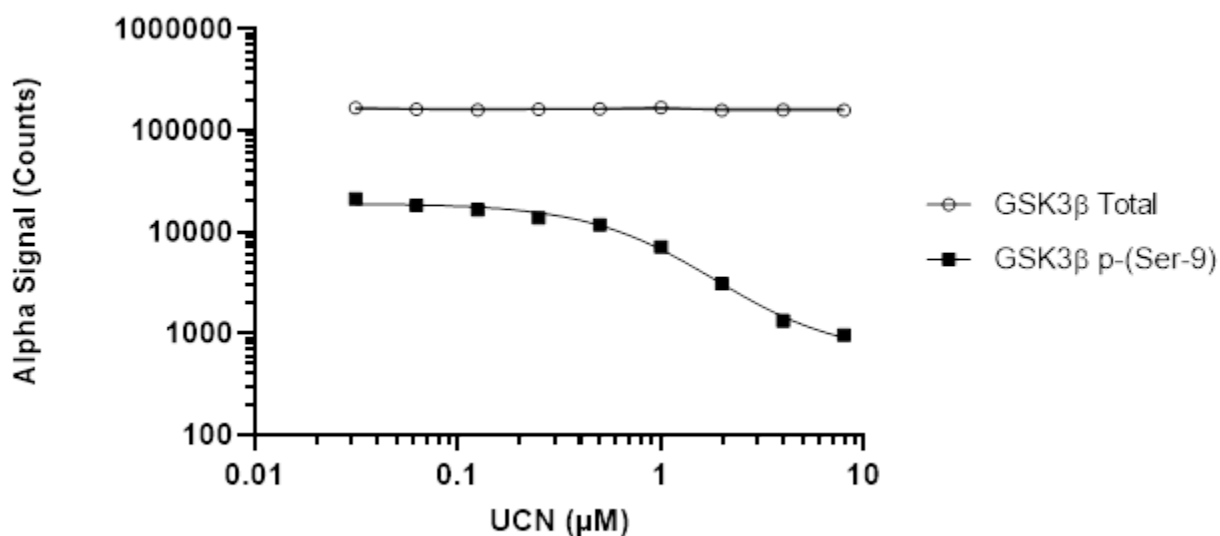
Positive Control Lysate: Prepared from MCF7 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 500 µg/mL Insulin for 30 minutes, lysed with 20 mL Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. SH-SY5Y cells were plated at 40K cells/well in a 96 well plate and incubated overnight. Cells were serum deprived for 2 hours then treated with Insulin at the indicated concentrations for 15 minutes. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Ser9) and Total GSK3β using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint.



Data obtained with a 2-plate, 2-incubation protocol. NIH3T3 cells were plated at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with UCN-01 at the indicated concentrations for 2 hours, then treated with 1.25 $\mu\text{g}/\text{mL}$ Insulin for 15 minutes. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Ser9) and Total GSK3 β using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint.



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, “Revvity”) do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product’s suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF”

