

AlphaLISA® SureFire® Ultra™ Terbium

Human and Mouse p-GSK3β (Ser9) Detection Kit

Product number: TBSU-PGSKB-A500, TBSU-PGSKB-A10K
TBSU-PGSKB-A50K, TBSU-PGSKB-A-HV



Kit specificity:

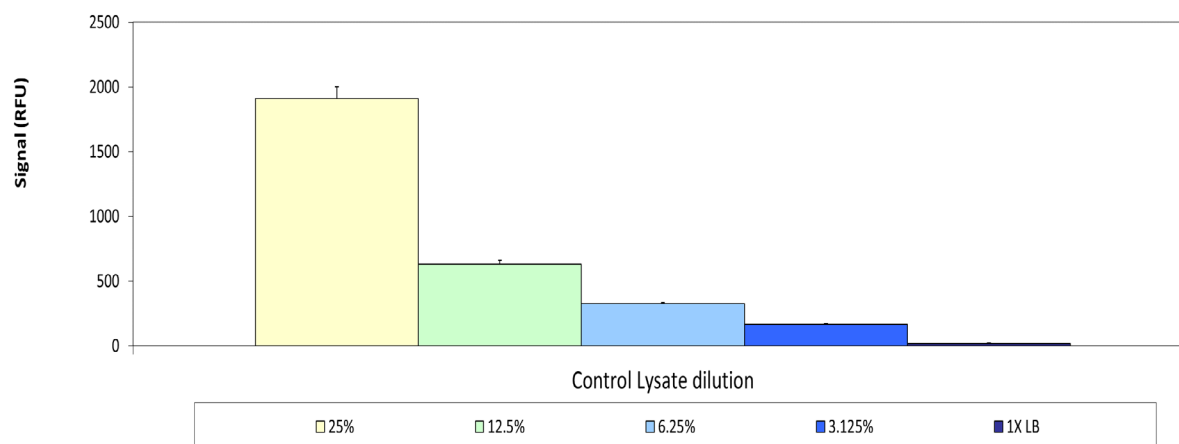
This assay kit contains antibodies which recognize the phospho-Ser9 epitope and a distal epitope on GSK3β. The protein detected by this kit corresponds to UniProt IDs P49841. GSK3β is also known as Glycogen synthase kinase-3 beta. These antibodies recognize GSK3β of human and mouse origin. Other species 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, treated with 500 µg/mL Insulin for 30 minutes and lysed with 20 mL of Lysis Buffer.

Representative data

Data obtained from the positive control lysate diluted in Lysis Buffer. Samples were assayed for Phospho GSK3β (Ser9) using the Terbium SureFire Ultra kit.



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"