

Anti-GSK3 α/β

Rabbit Polyclonal Antibody

Catalog # G09-63R-100

Lot # H245-8

Cited Applications

- Western blot (1:1000)

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the GSK3 α and GSK3 β protein

Cross Reactivity

- Western blot of GSK3 α and GSK3 β from human cells

GSK3 α and GSK3 β from other species may also be detectable

Host

Rabbit

Immunogen

Synthetic peptide corresponding to amino acids 214-218 of human GSK3 β

Formulation

PBS pH 7.4, 0.02% sodium azide, 50% glycerol

Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles.

Scientific Background

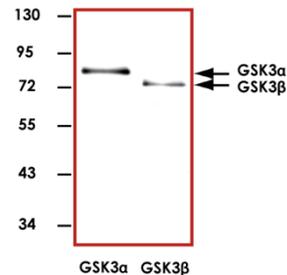
GSK3 α is a multifunctional protein serine kinase, homologous to Drosophila 'shaggy' (zeste-white3) and implicated in the control of several regulatory proteins including glycogen synthase and transcription factors (e.g., JUN) (1). GSK3 β is ubiquitously present in human tissues and implicated in the regulation of several physiological processes, including the control of glycogen and protein synthesis by insulin, modulation of the transcription factors AP-1 and CREB (2).

References

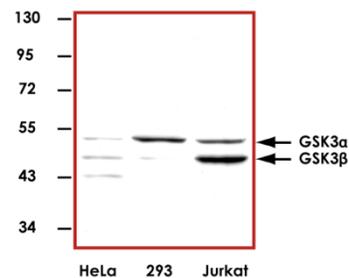
1. Ali, A. et al: Glycogen synthase kinase-3 : properties, functions, and regulation. Chem. Rev. 101: 2527-2540, 2001.
2. Sperber BR, Leight S, Goedert M, Lee VM. Glycogen synthase kinase-3 beta phosphorylates tau protein at multiple sites in intact cells. Neurosci Lett. 1995 Sep 8;197(2):149-53.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

Sample Data



Representative western blot with Anti-GSK3 α/β (1:1000) using 20 ng of GST tagged human recombinant GSK3 α and GSK3 β .



Representative western blot with Anti-GSK3 (1:1000) using 30 μ g of HeLa, HEK-293T, and Jurkat cell lysate.

Anti-GSK3 α/β

Rabbit Polyclonal Antibody

Catalog Number	G09-63R-100
Quantity	100 μ g
Specific Lot Number	H245-8
Purification	Affinity chromatography
Format	100 μ g in 100 μ l
Concentration	1.0 μ g/ μ l
Stability	1yr at -20°C from date of shipment
Storage & Shipping	Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.