
Phospho-GSK3B-Y216 Polyclonal Antibody

(Catalog #A72452)

Background

The protein encoded by this gene is a serine-threonine kinase, belonging to the glycogen synthase kinase subfamily. It is involved in energy metabolism, neuronal cell development, and body pattern formation. Polymorphisms in this gene have been implicated in modifying risk of Parkinson disease, and studies in mice show that overexpression of this gene may be relevant to the pathogenesis of Alzheimer disease. Alternatively spliced transcript variants encoding different isoforms have been found for this gene.

Description

Phospho-GSK3B-Y216 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.09% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P49841

Purification

Affinity Purification

Immunogen

A synthetic phosphorylated peptide around Y216 of human GSK3B (NP_001139628).

Storage

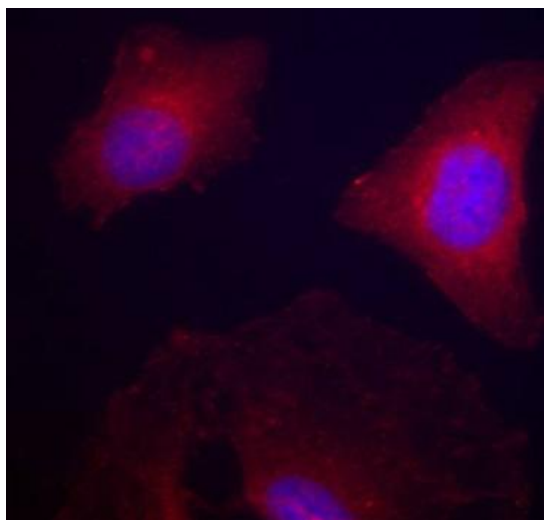
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

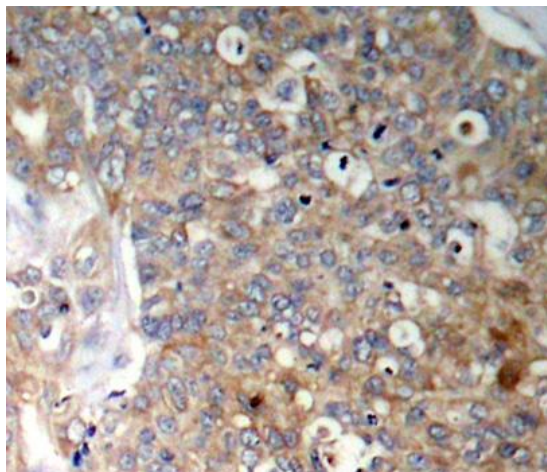
GSK3B; glycogen synthase kinase-3 beta

Application

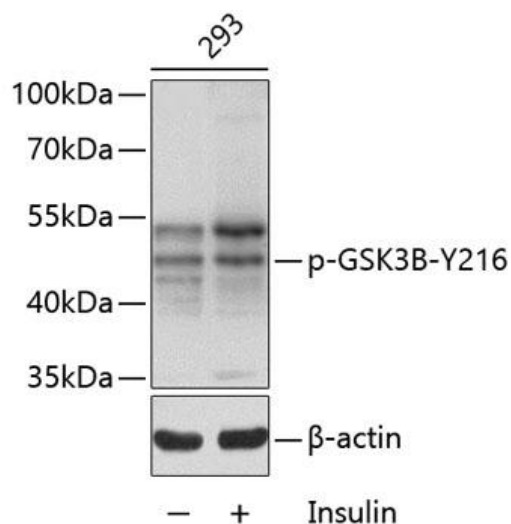
WB, IHC, IF, ELISA; Recommended dilution: WB 1:500 - 1:2000, IHC 1:50 - 1:100, IF 1:100 - 1:200, ELISA - Recommended starting concentration is 1 µg/mL. Please optimize the concentration based on your specific assay requirements.



Immunofluorescence analysis of methanol-fixed HeLa cells using Phospho-GSK3B-Y216 Polyclonal Antibody.



Immunohistochemistry of paraffin-embedded human breast carcinoma using Phospho-GSK3B-Y216 Polyclonal Antibody.



Western blot analysis of extracts of 293 cells, using Phospho-GSK3B-Y216 antibody at 1:1000 dilution. 293T cells were treated by Insulin (100nM) for 10 minutes after serum-starvation overnight. Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25ug per lane. Blocking buffer: 3% BSA.