[pSer21]GSK-3α monoclonal antibody (9B8)

Glycogen Synthase Kinase 3 (GSK3) exists in two isoforms, alpha and beta (GSK3 α , and GSK3 β). GSK3 is an active kinase that becomes phosphorylated by PKB/akt upon growth factor stimulation. Phosphorylation of GSK3 α at serine 21 or GSK3 β at serine 9, respectively, leads to inhibition of kinase activity of the GSK3.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

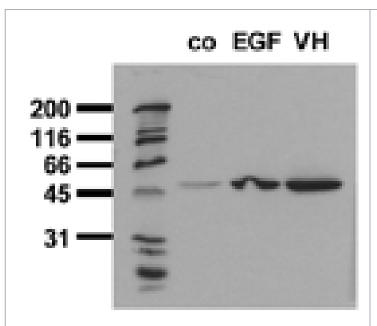
Order Online »

ADI-905-759-100

100µg

Manuals, SDS & CofA

View Online »



Western blot analysis of A431 cells untreated (co), EGF stimulated (EGF), or pervanadate treated (VH), with GSK3α (phospho-Ser21) mAb (9B8).

Handling & Storage

Use/Stability Reconstituted antibody is stable at -80°C up to 1 year, at 4°C up to 3 months.

Handling Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Glycogen synthate kinase-3α

Application ELISA, WB

Clone 9B8

Crossreactivity Does not cross-react with the homologous phosphorylation

site in GSK-3β (pSer9).

Formulation Lyophilized from 1ml of 2x PBS containing 0.09% sodium

azide, PEG, and sucrose.

Host Mouse

Immunogen Phosphopeptide conjugated to KLH.

Isotype IgG1

Purity Detail Thiophilic adsorption and size exclusion chromatography

purified.

Reconstitution Reconstitute with 1ml water (15 minutes at room

temperature).

Species Reactivity Dog, Human, Mouse, Rat

Specificity Recognizes GSK-3α phosphorylated at Ser²¹.

UniProt ID P49840

Last modified: May 29, 2024



info-

eu@enzolifesciences.com