

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

Glycogen Synthase Kinase 3 (GSK3) exists in two isoforms, alpha and beta (GSK3α, and GSK3β). GSK3 is an active kinase that becomes phosphorylated by PKB/akt upon growth factor stimulation. Phosphorylation of GSK3α at serine 21 or GSK3β at serine 9, respectively, leads to inhibition of kinase activity of the GSK3.

This antibody is covered by our [Worry-Free Guarantee](#).

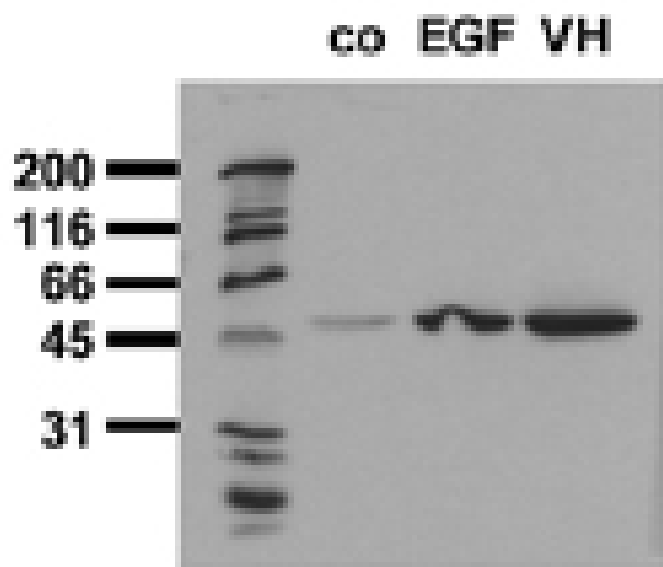
Ordering Information

[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 GSK3 α (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 Ser ²¹ .
UniProt ID	P49840

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)