

# Product Information



## GSK3 $\beta$ (Phospho-Ser<sup>9</sup>) Polyclonal Antibody

Item No. 10009374

<b>Synonym:</b>	Glycogen Synthase Kinase 3 $\beta$
<b>Supplied as:</b>	100 $\mu$ l of affinity-purified antibody in 10 mM HEPES (pH 7.5), 150 mM NaCl, 100 $\mu$ g BSA per ml and 50% glycerol.
<b>Host:</b>	Rabbit
<b>Antigen:</b>	Phosphopeptide corresponding to amino acid residues surrounding the phospho-Ser <sup>9</sup> of GSK3 $\beta$
<b>Cross Reactivity:</b>	(+) Rat GSK3 $\beta$ ; expected to react with human, mouse, chicken, bovine, canine, non-human primate, zebrafish, and <i>Xenopus</i> GSK3 $\beta$
<b>Stability:</b>	$\geq$ 1 year at -20°C
<b>Application:</b>	The recommended starting dilution for western blot is 1:1,000. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

Glycogen synthase kinase 3 (GSK3) is a serine/threonine kinase that is involved in the regulation of many signaling pathways. To date, two isoforms have been identified: GSK3 $\alpha$  and GSK3 $\beta$ . Specifically, GSK3 $\beta$  has been shown to play a key inhibitory role in both the insulin and Wnt signaling pathways.<sup>1</sup> It has been suggested that Ser<sup>9</sup> phosphorylation underlies the inhibition of GSK3 $\beta$  by insulin.<sup>2</sup>

### References

1. Papkoff, J. and Aikawa, M. WNT-1 and HGF regulate GSK3 $\beta$  activity and  $\beta$  catenin signaling in mammary epithelial cells. *Biochem. Biophys. Res. Commun.* **247(3)**, 851-858 (1998).
2. Sutherland, C., Leighton, I.A., and Cohen, P. Inactivation of glycogen synthase kinase 3  $\beta$  by phosphorylation: New kinase connections in insulin and growth factor signaling. *Biochem J.* **296(Pt. 1)**, 15 (1993).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/10009374](http://www.caymanchem.com/catalog/10009374)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 12/20/2011

### Cayman Chemical

#### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

#### Phone

(800) 364-9897  
(734) 971-3335

#### Fax

(734) 971-3640

#### E-Mail

[custserv@caymanchem.com](mailto:custserv@caymanchem.com)

#### Web

[www.caymanchem.com](http://www.caymanchem.com)