

PRODUCT INFORMATION



GSK3β Inhibitor VIII

Item No. 14704

CAS Registry No.: 487021-52-3

Formal Name: N-[(4-methoxyphenyl)methyl]-N'-(5-nitro-2-thiazolyl)-urea

Synonyms: AR 014418, AR 0133418, Glycogen Synthase Kinase 3β Inhibitor VIII, SN 4521

MF: C₁₂H₁₂N₄O₄S

FW: 308.3

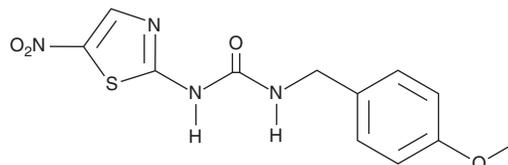
Purity: ≥98%

UV/Vis.: λ_{max}: 226, 275, 348, 353 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

GSK3β inhibitor VIII is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK3β inhibitor VIII in the solvent of choice, which should be purged with an inert gas. GSK3β inhibitor VIII is soluble in the organic solvent DMSO at a concentration of approximately 30 mg/ml.

Description

GSK3β inhibitor VIII is an ATP-competitive inhibitor of the serine/threonine kinase, glycogen synthase kinase 3β ($K_i = 38$ nM; $IC_{50} = 104$ nM).¹ It does not significantly inhibit the closely related cyclin-dependent kinase (Cdk)2 or Cdk5 ($K_{iS} = > 100$ μM).¹ GSK3β inhibitor VIII demonstrates neuroprotective effects by inhibiting tau phosphorylation ($IC_{50} = 2.7$ μM) at a GSK3-specific site (Ser³⁹⁶) in cells stably expressing human four-repeat tau protein and also by protecting β-amyloid-exposed N2A neuroblastoma cells against cell death.¹

Reference

1. Bhat, R., Xue, Y., Berg, S., *et al.* Structural insights and biological effects of glycogen synthase kinase 3-specific inhibitor AR-A014418. *J. Biol. Chem.* **278**(46), 45937-45945 (2003).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/23/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM