

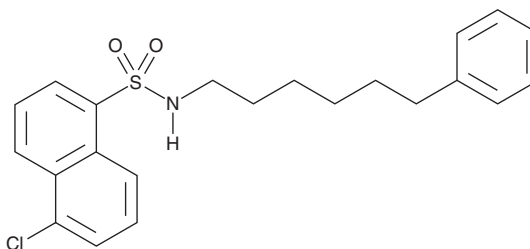
PRODUCT INFORMATION



SC-9

Item No. 30910

CAS Registry No.: 102649-78-5
Formal Name: 5-chloro-N-(6-phenylhexyl)-1-naphthalenesulfonamide
MF: C₂₂H₂₄ClNO₂S
FW: 402.0
Purity: ≥95%
Supplied as: A 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

SC-9 is supplied as a solid. A stock solution may be made by dissolving the SC-9 in the solvent of choice, which should be purged with an inert gas. SC-9 is slightly soluble in acetonitrile and chloroform.

Description

SC-9 is an activator of protein kinase C (PKC).¹ It is selective for PKC over myosin light-chain kinase (MLCK), PKA, and PKG at 300 μM. SC-9 (200 μM) increases PKC-induced MLC phosphorylation in the presence of calcium.³ It induces proliferation of quiescent primary neonatal mouse epidermal melanoblasts when used at a concentration of 10 μM.² SC-9 (2.5-20 μM) reduces the motility of isolated intact chicken sperm.⁴

References

1. Ito, M., Tanaka, T., Inagaki, M., *et al.* N-(6-phenylhexyl)-5-chloro-1-naphthalenesulfonamide, a novel activator of protein kinase C. *Biochemistry* **25(15)**, 4179-4184 (1986).
2. Hirobe, T. Effects of activators (SC-9 and OAG) and inhibitors (staurosporine and H-7) of protein kinase C on the proliferation of mouse epidermal melanoblasts in serum-free culture. *J. Cell Sci.* **107(Pt 6)**, 1679-1686 (1994).
3. Nishino, H., Kitagawa, K., Iwashima, A., *et al.* N-(6-Phenylhexyl)-5-chloro-1-naphthalenesulfonamide is one of a new class of activators for Ca²⁺-activated, phospholipid-dependent protein kinase. *Biochim. Biophys. Acta* **889(2)**, 236-239 (1986).
4. Ashizawa, K., Katayama, S., Kobayashi, T., *et al.* Possible role of protein kinase C in regulation of flagellar motility and intracellular free Ca²⁺ concentration of fowl spermatozoa. *J. Reprod. Fertil.* **101(3)**, 511-517 (1994).

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, 10/20/2022

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