

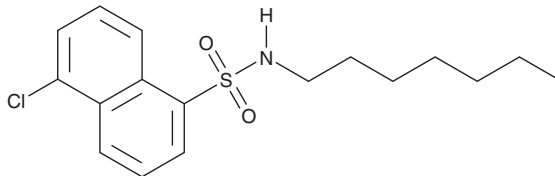
PRODUCT INFORMATION



SC-10

Item No. 42947

CAS Registry No.: 102649-79-6
Formal Name: 5-chloro-N-heptyl-1-naphthalenesulfonamide
MF: $C_{17}H_{22}ClNO_2S$
FW: 339.9
Purity: $\geq 98\%$
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

SC-10 is supplied as a crystalline solid. A stock solution may be made by dissolving the SC-10 in the solvent of choice, which should be purged with an inert gas. SC-10 is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

SC-10 is an activator of PKC.¹ It induces histone phosphorylation by recombinant human PKC in a cell-free assay when used at concentrations ranging from 100 to 500 μM .

Reference

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).

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, 03/13/2025

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