

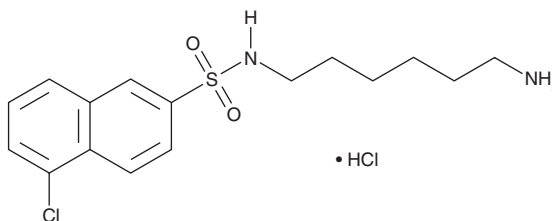
PRODUCT INFORMATION



W-9 (hydrochloride)

Item No. 43152

CAS Registry No.: 69762-85-2
Formal Name: N-(6-aminohexyl)-5-chloro-2-naphthalenesulfonamide, monohydrochloride
MF: $C_{16}H_{21}ClN_2O_2S \cdot HCl$
FW: 377.3
Purity: $\geq 98\%$
Supplied as: A 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

W-9 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the W-9 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. W-9 (hydrochloride) is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

W-9 is a calmodulin antagonist ($IC_{50} = 16 \mu M$ in a radioligand binding assay).¹ It inhibits calmodulin-dependent phosphodiesterase (PDE) activation in a cell-free assay ($IC_{50} = 14 \mu M$) and erythrocyte ghost ATPase activation ($IC_{50} = 15-18 \mu M$).^{1,2} W-9 induces relaxation in isolated rabbit aortic strips precontracted with prostaglandin $F_{2\alpha}$ ($PGF_{2\alpha}$; $EC_{50} = 72 \mu M$) and inhibits ATP-induced superprecipitation of bovine aorta smooth muscle actomyosin ($IC_{50} = 46 \mu M$).³ It also inhibits the replication of dengue virus in HEK293 cells ($IC_{50} = 3.19 \mu M$).⁴

References

1. Hidaka, H. and Tanaka, T. Naphthalenesulfonamides as calmodulin antagonists. *Methods Enzymol.* **102**, 185-194 (1983).
2. Kobayashi, R., Tawata, M., and Hidaka, H. Ca^{2+} regulated modulator protein interacting agents: inhibition of Ca^{2+} - Mg^{2+} -ATPase of human erythrocyte ghost. *Biochem. Biophys. Res. Commun.* **88(3)**, 1037-1045 (1979).
3. Hidaka, H., Yamaki, T., Totsuka, T., *et al.* Selective inhibitors of Ca^{2+} -binding modulator of phosphodiesterase produce vascular relaxation and inhibit actin-myosin interaction. *Mol. Pharmacol.* **15(1)**, 49-59 (1979).
4. Shum, D., Smith, J.L., Hirsch, A.J., *et al.* High-content assay to identify inhibitors of dengue virus infection. *Assay Drug Dev. Technol.* **8(5)**, 553-570 (2010).

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, 04/15/2025

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