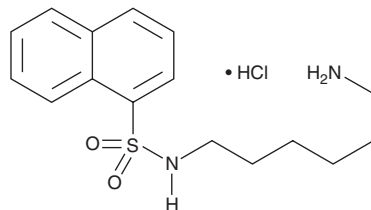


PRODUCT INFORMATION

W-5 (hydrochloride)

Item No. 14271

CAS Registry No.: 61714-25-8
Formal Name: N-(6-aminohexyl)-1-naphthalenesulfonamide, monohydrochloride
MF: C₁₆H₂₂N₂O₂S • HCl
FW: 342.9
Purity: ≥98%
UV/Vis.: λ_{max}: 223, 289 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

W-5 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the W-5 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. W-5 (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of W-5 (hydrochloride) in ethanol is approximately 0.33 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of W-5 (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of W-5 (hydrochloride) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

W-5 is a chlorine-deficient analog of W-7 (Item No. 14826), a calmodulin antagonist that binds to calmodulin and inhibits Ca²⁺/calmodulin-regulated enzyme activities (K_i = 11 μM).^{1,2} W-5 is less active than its calmodulin antagonist counterpart, inhibiting CHO-K1, MCF-7 breast cancer, and human myeloid progenitor cell proliferation with IC₅₀ values greater than 100 μM.^{1,3,4} As a weaker antagonist for calmodulin, W-5 is suitable for use as a control compound for understanding the specificity of other calmodulin antagonists.

References

1. Hidaka, H., Sasaki, Y., Tanaka, T., *et al.* N-(6-aminoethyl)-5-chloro-1-naphthalenesulfonamide, a calmodulin antagonist, inhibits cell proliferation. *Proc. Natl. Acad. Sci. USA* **78**(7), 4354-4357 (1981).
2. Tanaka, T. and Hidaka, H. Hydrophobic regions function in calmodulin-enzyme(s) interactions. *J. Biol. Chem.* **255**(23), 11078-11080 (1980).
3. Gulino, A., Barrera, G., Vacca, A., *et al.* Calmodulin antagonism and growth-inhibiting activity of triphenylethylene antiestrogens in MCF-7 human breast cancer cells. *Cancer Res.* **46**, 6274-6278 (1986).
4. Katayama, N., Nishikawa, M., Komada, F., *et al.* A role for calmodulin in the growth of human hematopoietic progenitor cells. *Blood* **75**(7), 1446-1454 (1990).

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, 10/21/2022

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