

PRODUCT INFORMATION



Noscapine (hydrochloride)

Item No. 17255

CAS Registry No.: 912-60-7
Formal Name: (3S)-6,7-dimethoxy-3-[(5R)-5,6,7,8-tetrahydro-4-methoxy-6-methyl-1,3-dioxolo[4,5-g]isoquinolin-5-yl]-1(3H)-isobenzofuranone, monohydrochloride

Synonyms: Narcotine, Opianine

MF: C₂₂H₂₃NO₇ • HCl

FW: 449.9

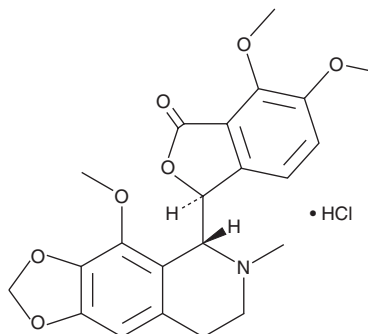
Purity: ≥98%

UV/Vis.: λ_{max}: 214, 313 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.

Description

Noscapine is an opioid alkaloid with antitussive and antimitotic properties.¹ It binds to tubulin and arrests mammalian cells in mitosis at a concentration of 20 μM.² Noscapine induces apoptosis (IC₅₀ = 25 and 10 μM in HeLa cells and thymocytes, respectively) and demonstrates antitumor activity both *in vitro* (IC₅₀ = 42.3 and 39.1 μM in MCF-7 and Renal 1983 cell lines, respectively) and in mice.²

References

1. Kingston, D.G.I. Tubulin-interactive natural products as anticancer agents. *J. Nat. Prod.* **72**(3), 507-515 (2009).
2. Ye, K., Ke, Y., Keshava, N., *et al.* Opium alkaloid noscapine is an antitumor agent that arrests metaphase and induces apoptosis in dividing cells. *Proc. Natl. Acad. Sci. USA* **95**(4), 1601-1606 (1998).

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, 01/03/2023

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