

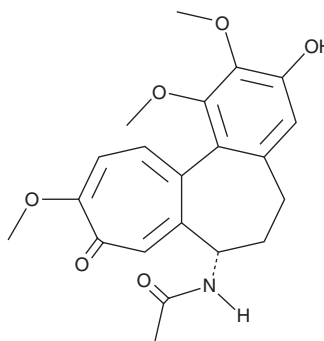
PRODUCT INFORMATION



3-demethyl Colchicine

Item No. 35238

CAS Registry No.: 7336-33-6
Formal Name: N-[(7S)-5,6,7,9-tetrahydro-3-hydroxy-1,2,10-trimethoxy-9-oxobenzo[a]heptalen-7-yl]-acetamide
Synonym: (-)-3-Demethylcolchicine
MF: C₂₁H₂₃NO₆
FW: 385.4
Purity: ≥98%
UV/Vis.: λ_{max}: 242, 352 nm
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

3-demethyl Colchicine is supplied as a solid. A stock solution may be made by dissolving the 3-demethyl colchicine in the solvent of choice, which should be purged with an inert gas. 3-demethyl Colchicine is soluble in ethanol.

Description

3-demethyl Colchicine is an active metabolite of colchicine (Item No. 9000760).^{1,2} It suppresses antibody production induced by sheep red blood cells in co-culture with isolated mouse spleen cells when used at concentrations ranging from 0.1 to 10 µg/ml.² 3-demethyl Colchicine (100 µg/animal) inhibits carrageenan-induced paw edema in rats.³

References

- Schönharting, M., Mende, G., and Siebert, G. Metabolic transformation of colchicine. II. The metabolism of colchicine by mammalian liver microsomes. *Hoppe Seylers Z. Physiol. Chem.* **355(11)**, 1391-1399 (1974).
- Sterzl, J., Santavý, F., Sedmera, P., et al. Effect of colchicine derivatives on the antibody response induced in vitro. *Folia Microbiol. (Praha)* **27(4)**, 256-266 (1982).
- Sugio, K., Maruyama, M., Tsurufuji, S., et al. Separation of tubulin-binding and anti-inflammatory activity in colchicine analogs and congeners. *Life Sci.* **40(1)**, 35-39 (1987).

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, 12/12/2022

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