

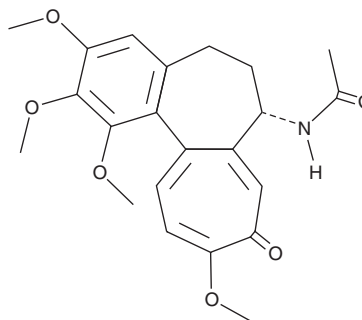
PRODUCT INFORMATION



Colchicine

Item No. 9000760

CAS Registry No.: 64-86-8
Formal Name: N-[(7S)-5,6,7,9-tetrahydro-1,2,3,10-tetramethoxy-9-oxobenzo[a]heptalen-7-yl]-acetamide
MF: C₂₂H₂₅NO₆
FW: 399.4
Purity: ≥98%
UV/Vis.: λ_{max}: 202, 243, 352 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

Colchicine is supplied as a crystalline solid. A stock solution may be made by dissolving the colchicine in the solvent of choice, which should be purged with an inert gas. Colchicine is soluble in organic solvents such as ethanol and DMSO. The solubility of colchicine in these solvents is approximately 25 mg/ml.

Description

Colchicine is an inhibitor of microtubule polymerization (IC₅₀ = 3.2 μM) that binds to tubulin, which disrupts spindle formation during mitosis.¹ It inhibits growth of MCF-7 human breast carcinoma cells with an IC₅₀ value of 13 nM.¹ Colchicine has anti-inflammatory activity, inhibiting neutrophil motility and activity when used at a dose of 5 μmol/kg in a mouse model of gout and preventing the deposition of uric acid.^{2,3}

References

1. Martino, G.D., Regina, G.L., Coluccia, A., *et al.* Arylthioindoles, potent inhibitors of tubulin polymerization. *J. Med. Chem.* **47(25)**, 6120-6123 (2004).
2. McCarty, D.J. Urate crystals, inflammation, and colchicine. *Arthritis Rheum.* **58(2)**, S20-S24 (2008).
3. Chia, E.W., Grainger, R., and Harper, J.L. Colchicine suppresses neutrophil superoxide production in a murine model of gouty arthritis: A rationale for use of low-dose colchicine. *Br. J. Pharmacol.* **153(6)**, 1288-95 (2008).

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, 11/08/2022

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