

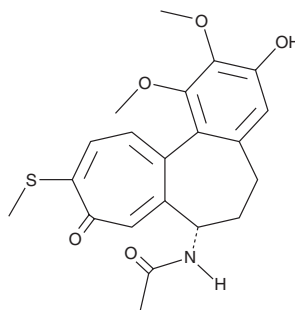
PRODUCT INFORMATION



3-O-demethyl Thiocolchicine

Item No. 35497

CAS Registry No.: 87424-25-7
Formal Name: N-[(7S)-5,6,7,9-tetrahydro-3-hydroxy-1,2-dimethoxy-10-(methylthio)-9-oxobenzo[a]heptalen-7-yl]-acetamide
Synonyms: NSC 361792, 3-Demethylthiocolchicine
MF: $C_{21}H_{23}NO_5S$
FW: 401.5
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 257, 386 nm
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

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

Description

3-O-demethyl Thiocolchicine is a degradation product of the spasmolytic agent thiocolchicoside.¹ It is formed from thiocolchicoside under acidic and basic conditions.

Reference

1. Erika, D.G., Silvio, A., and Giorgio, G. Forced degradation study of thiocolchicoside: Characterization of its degradation products. *J. Pharm. Biomed. Anal.* **61**, 215-223 (2012).

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