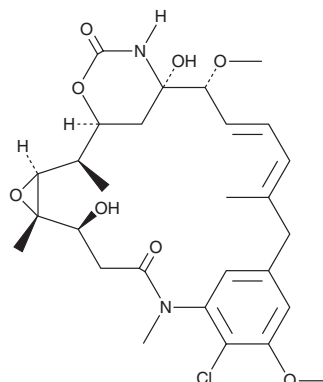


# PRODUCT INFORMATION



## Maytansinol Item No. 29228

**CAS Registry No.:** 57103-68-1  
**Formal Name:** 3-O-de[2-(acetylmethylamino)-1-oxopropyl]-maytansine  
**Synonyms:** Ansamitocin P-O, Antibiotic C-15003P0, NSC 239386  
**MF:** C<sub>28</sub>H<sub>37</sub>ClN<sub>2</sub>O<sub>8</sub>  
**FW:** 565.1  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 232, 244, 253 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Synthetic



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Maytansinol is supplied as a crystalline solid. A stock solution may be made by dissolving the maytansinol in the solvent of choice, which should be purged with an inert gas. Maytansinol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of maytansinol in these solvents is approximately 10, 20, and 25 mg/ml, respectively.

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 maytansinol can be prepared by directly dissolving the crystalline solid in aqueous buffers. Maytansinol is slightly soluble in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

### Description

Maytansinol is an ansa macrolide originally isolated from *P. verrucosa* that has antimitotic and anticancer activities.<sup>1-3</sup> It inhibits polymerization and induces depolymerization of bovine brain tubulin with EC<sub>50</sub> values of 12 and 43 μM, respectively.<sup>2</sup> Maytansinol inhibits sea urchin egg mitosis when used at a concentration of 10 μM and decreases proliferation of KB nasopharyngeal cancer cells (EC<sub>50</sub> = 0.19 μg/ml).<sup>3</sup>

### References

1. Kupchan, S.M., Branfman, A.R., Sneden, A.T., *et al.* Novel maytansinoids. Naturally occurring and synthetic antileukemic esters of maytansinol. *J. Am. Chem. Soc.* **97(18)**, 5294-5295 (1975).
2. Ikeyama, S. and Takeuchi, M. Antitubulin activities of ansamitocins and maytansinoids. *Biochem. Pharmacol.* **30(17)**, 2421-2425 (1981).
3. Kupchan, S.M., Sneden, A.T., Branfman, A.R., *et al.* Structural requirements for antileukemic activity among the naturally occurring and semisynthetic maytansinoids. *J. Med. Chem.* **21(1)**, 31-37 (1978).

#### 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, 12/13/2022

#### 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