

PRODUCT INFORMATION



Val-Cit-PAB-MMAD

Item No. 45988

CAS Registry No.: 1415329-13-3

Formal Name: N-methyl-N-[[[4-[[L-valyl-N-(aminocarbonyl)-L-ornithyl]amino]phenyl]methoxy]carbonyl]-L-valyl-N-[(1S,2R)-2-methoxy-4-[(2S)-2-[(1R,2R)-1-methoxy-2-methyl-3-oxo-3-[[[(1S)-2-phenyl-1-(2-thiazolyl)ethyl]amino]propyl]-1-pyrrolidinyl]-1-[(1S)-1-methylpropyl]-4-oxobutyl]-N-methyl-L-valinamide

Synonyms: Val-Cit-PAB-monomethyl Auristatin D, Val-Cit-PAB-monomethyl Dolastatin 10

MF: $C_{60}H_{93}N_{11}O_{11}S$

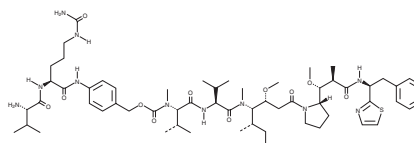
FW: 1,176.5

Purity: $\geq 98\%$

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

Val-Cit-PAB-MMAD is supplied as a solid. A stock solution may be made by dissolving the Val-Cit-PAB-MMAD in the solvent of choice, which should be purged with an inert gas. Val-Cit-PAB-MMAD is sparingly soluble (1-10 mg/ml) in a 1:1 solution of acetonitrile:water.

Description

Val-Cit-PAB-MMAD is a building block that contains the tubulin inhibitor monomethyl auristatin D (MMAD) and a Val-Cit-PAB linker region. It has been used in the synthesis of antibody-drug conjugates (ADCs) containing trastuzumab that are cytotoxic to breast cancer cells *in vitro*.¹

Reference

1. Chen, B., Gianolio, D.A., Stefano, J.E., et al. Design, synthesis, and in vitro evaluation of multivalent drug linkers for high-drug-load antibody-drug conjugates. *ChemMedChem* **13**(8), 790-794 (2018).

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, 07/01/2026

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