

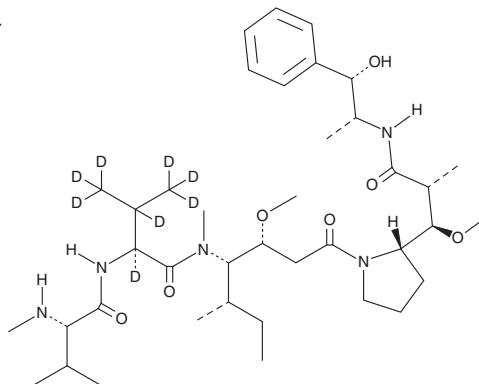
PRODUCT INFORMATION



Monomethyl Auristatin E-d₈

Item No. 44386

CAS Registry No.: 2070009-72-0
Formal Name: N-methyl-L-valyl-N-[(1S,2R)-4-[(2S)-2-[(1R,2R)-3-[[[(1R,2S)-2-hydroxy-1-methyl-2-phenylethyl]amino]-1-methoxy-2-methyl-3-oxopropyl]-1-pyrrolidinyl]-2-methoxy-1-[(1S)-1-methylpropyl]-4-oxobutyl]-N-methyl-L-valinamide-2,3,4,4,4,5,5,5-d₈
Synonyms: Brentuximab Vedotin-d₈, MMAE-d₈
MF: C₃₉H₅₉D₈N₅O₇
FW: 726.0
Chemical Purity: ≥95% (Monomethyl Auristatin E)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₈); ≤1% d₀
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

Monomethyl auristatin E-d₈ is intended for use as an internal standard for the quantification of monomethyl auristatin E (MMAE; Item No. 16267) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

MMAE-d₈ is supplied as a solid. A stock solution may be made by dissolving the MMAE-d₈ in the solvent of choice, which should be purged with an inert gas. MMAE-d₈ is soluble in acetonitrile, methanol, DMSO, and dimethyl formamide.

Description

MMAE is a derivative of dolastatin 10 and an inhibitor of tubulin polymerization.¹ It inhibits the assembly of MAP-rich tubulin and alters the morphology of microtubule ends in cell-free assays when used at concentrations ranging from 1 to 10 μM. MMAE induces cell cycle arrest at the G₂/M phase in, and inhibits the proliferation of, MCF-7 cells (IC₅₀s = 0.5 and 0.9 nM, respectively). It is cytotoxic against a panel of nine lymphoma cell lines (IC₅₀s = 0.039-1.35 nM).² MMAE has been used as the cytotoxic payload component of antibody-drug conjugates (ADCs) with anticancer activity *in vitro* and in animal models.^{1,2}

References

1. Best, R.L., LaPointe, N.E., Azarenko, O., *et al.* Microtubule and tubulin binding and regulation of microtubule dynamics by the antibody drug conjugate (ADC) payload, monomethyl auristatin E (MMAE): Mechanistic insights into MMAE ADC peripheral neuropathy. *Toxicol. Appl. Pharmacol.* **421**, 115534 (2021).
2. Francisco, J.A., Cervený, C.G., Meyer, D.L., *et al.* cAC10-vcMMAE, an anti-CD30-monomethyl auristatin E conjugate with potent and selective antitumor activity. *Blood* **102(4)**, 1458-1465 (2003).

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, 09/04/2025

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