

PRODUCT INFORMATION



VcMMAE

Item No. 35646

CAS Registry No.: 646502-53-6

Formal Name: N-[[[4-[[N-[6-(2,5-dihydro-2,5-dioxo-1H-pyrrol-1-yl)-1-oxohexyl]-L-valyl-N⁵-(aminocarbonyl)-L-ornithyl]amino]phenyl]methoxy]carbonyl]-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-pyrrolidiny]-2-methoxy-1-[(1S)-1-methylpropyl]-4-oxobutyl]-N-methyl-L-valinamide

Synonyms: MC-VC-PAB-MMAE, Vedotin

MF: C₆₈H₁₀₅N₁₁O₁₅

FW: 1,316.6

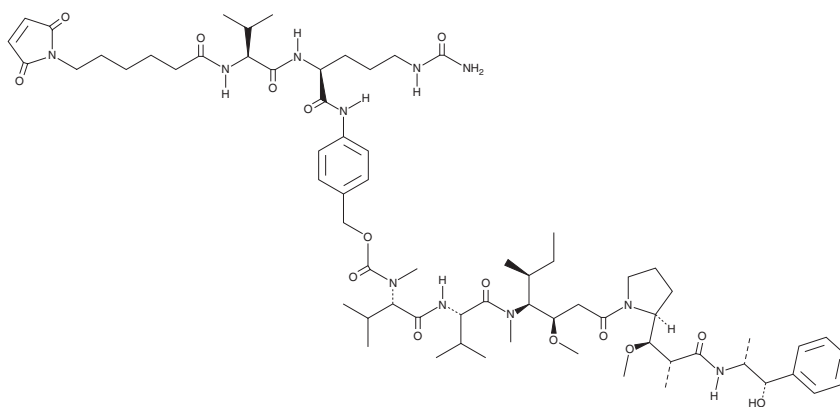
Purity: ≥98%

UV/Vis.: λ_{max}: 251 nm

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

VcMMAE is supplied as a solid. A stock solution may be made by dissolving the VcMMAE in the solvent of choice, which should be purged with an inert gas. VcMMAE is soluble in ethanol and DMSO.

Description

VcMMAE is a drug-linker molecule that has been used in the synthesis of antibody-drug conjugates (ADCs).^{1,2} It contains a maleimidocaproyl (Mc) group that can be conjugated to an antibody, a protease-cleavable peptide linker, and the antimitotic and anticancer agent monomethyl auristatin E (MMAE; Item No. 16267).

References

1. Chuprakov, S., Ogunkoya, A.O., Barfield, R.M., *et al.* Tandem-cleavage linkers improve the *in vivo* stability and tolerability of antibody-drug conjugates. *Bioconj. Chem.* **32**(4), 746-754 (2021).
2. Dimou, M., Papageorgiou, S.G.S., N., Katodritou, E., *et al.* Real-life experience with the combination of polatuzumab vedotin, rituximab, and bendamustine in aggressive B-cell lymphomas. *Hematol. Oncol.* **39**(3), 336-348 (2021).

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