

PRODUCT INFORMATION



Val-Cit-PAB-MMAE

Item No. 41048

CAS Registry No.: 644981-35-1

Formal Name: N-methyl-N-[[[4-[[L-valyl-N⁵-(aminocarbonyl)-L-ornithyl]amino]phenyl]methoxy]carbonyl]-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

Synonyms: N-PM-0008, VC-PAB-MMAE

MF: C₅₈H₉₄N₁₀O₁₂

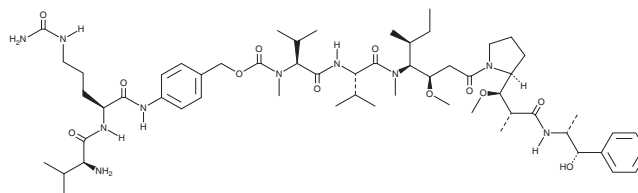
FW: 1,123.4

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

Description

Val-Cit-PAB-MMAE is a building block that has been used in the synthesis of drug-linker molecules for antibody-drug conjugates (ADCs).¹⁻³

References

1. 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).
2. Tse, K.F., Jeffers, M., Pollack, V.A., *et al.* CR011, a fully human monoclonal antibody-auristatin E conjugate for the treatment of melanoma. *Clin. Cancer Res.* **12**(4), 1373-1382 (2006).
3. Chuprakov, S., Ogunkoya, A.O., Barfield, R.M., *et al.* Tandem-cleavage linkers improve the *in vivo* stability and tolerability of antibody-drug conjugates. *Bioconjug. Chem.* **32**(4), 746-754 (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, 05/31/2024

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