

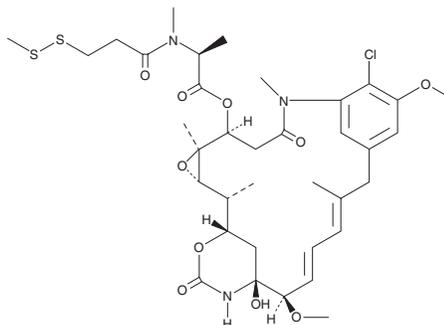
PRODUCT INFORMATION



DM1-SMe

Item No. 22477

CAS Registry No.: 138148-68-2
Formal Name: N²-deacetyl-N²-[3-(methylthio)-1-oxopropyl]-maytansine
MF: C₃₆H₅₀ClN₃O₁₀S₂
FW: 784.4
Purity: ≥98%
UV/Vis.: λ_{max}: 234, 254, 280 nm
Supplied as: A crystalline 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

DM1-SMe is supplied as a crystalline solid. A stock solution may be made by dissolving the DM1-SMe in the solvent of choice, which should be purged with an inert gas. DM1-SMe is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of DM1-SMe in DMSO is approximately 12 mg/ml and approximately 16 mg/ml in DMF.

DM1-SMe is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, DM1-SMe should first be dissolved in DMF and then diluted with the aqueous buffer of choice. DM1-SMe has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

DM1-SMe is an analog of mertansine (DM1; Item No. 22483) with a thiomethane cap attached to the sulfhydryl group.¹ It has *in vitro* activity against the pediatric preclinical testing program (PPTP) panel of cancer cell lines with IC₅₀ values of 0.002 to >3 nM.

References

1. Wood, A.C., Maris, J.M., Gorlick, R., *et al.* Initial testing (Stage 1) of the antibody-maytansinoid conjugate, IMG901 (Lorvotuzumab mertansine), by the pediatric preclinical testing program. *Pediatr. Blood Cancer* **60(11)**, 1860-1867 (2013).

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, 11/04/2022

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