

PRODUCT INFORMATION



MMAF

Item No. 30197

CAS Registry No.: 745017-94-1
Formal Name: N-methyl-L-valyl-L-valyl-(3R,4S,5S)-3-methoxy-5-methyl-4-(methylamino)heptanoyl-(α R, β R,2S)- β -methoxy- α -methyl-2-pyrrolidinepropanoyl-L-phenylalanine

Synonym: Monomethylauristatin F

MF: C₃₉H₆₅N₅O₈

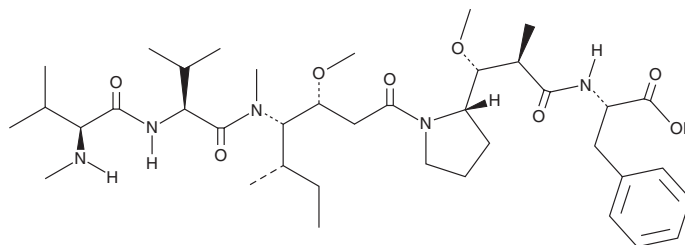
FW: 732.0

Purity: \geq 95%

Supplied as: A solid

Storage: -20°C

Stability: \geq 2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

MMAF is supplied as a solid. A stock solution may be made by dissolving the MMAF in the solvent of choice, which should be purged with an inert gas. MMAF is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of MMAF in these solvents is approximately 2, 25, and 10 mg/ml, respectively.

MMAF is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MMAF should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. MMAF has a solubility of approximately 0.14 mg/ml in a 1:6 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

MMAF is an auristatin with anticancer activity.¹ It is cytotoxic to Karpas 299 anaplastic large cell lymphoma, H3396 breast carcinoma, and 786-O and Caki-1 renal carcinoma cells *in vitro* (IC₅₀s = 119, 105, 257, and 200 nM, respectively), as well as in a panel of 11 lymphoma cell lines (IC₅₀s = 34.8-164 nM).

Reference

1. Doronina, S.O., Mendelsohn, B.A., Bovee, T.D., *et al.* Enhanced activity of monomethylauristatin F through monoclonal antibody delivery: Effects of linker technology on efficacy and toxicity. *Biconjug. Chem.* **17**(1), 114-124 (2006).

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, 05/21/2020

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