

PRODUCT INFORMATION



MMAE Monoclonal Antibody (Clone 1G1)

Item No. 44706

Overview and Properties

Contents:	This vial contains 100 µg of protein G-purified monoclonal antibody.
Synonyms:	Monomethyl Auristatin E, Vedotin
Immunogen:	Vc-MMAE-KLH
Cross Reactivity:	(+++ MMAE, (+) Mertansine
Species Reactivity:	(+) Species independent
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	PBS (pH 7.2), with 50% glycerol and 0.02% sodium azide
Clone:	1G1
Host:	Mouse
Isotype:	IgG _{3k}
Applications:	ELISA and Western blot (WB); the recommended starting dilutions are 1:1,000 (2 µg/ml) for both WB and ELISA. Other applications were not tested; therefore, optimal working concentration/dilution should be determined empirically.

Image

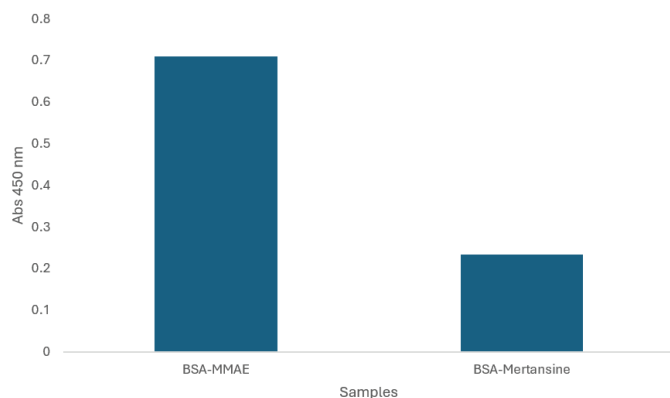


Figure 1. MMAE Monoclonal Antibody (Clone 1G1) (Item No. 44706) detected an MMAE-albumin conjugate with 32% cross-reactivity to a mertansine-albumin conjugate by ELISA.

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, 01/15/2026

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

PRODUCT INFORMATION



Description

Monomethyl auristatin E (MMAE) is an inhibitor of tubulin polymerization and a derivative of dolastatin 10.¹ It inhibits the assembly of MAP-rich tubulin and induces cell cycle arrest and cell death in cancer cells *in vitro*.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} Cayman's MMAE Monoclonal Antibody (Clone 1G1) can be used for ELISA and Western blot (WB) applications.

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).

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