

PRODUCT INFORMATION



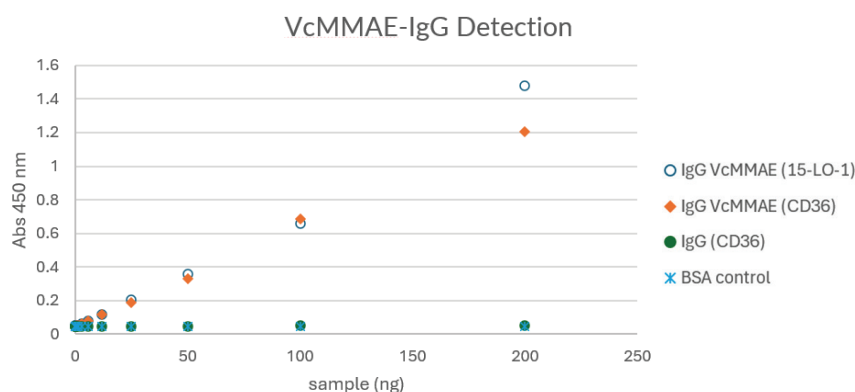
VcMMAE Polyclonal Antibody - Biotinylated

Item No. 42295

Overview and Properties

Contents:	This vial contains 200 µg of protein A-affinity purified polyclonal antibody.
Synonyms:	VC-MMAE, VC-Monomethyl Auristatin E Conjugate
Immunogen:	VcMMAE-KLH conjugate
Cross Reactivity:	(+) VcMMAE-labeled antibodies; (-) Albumin, Unlabeled antibodies
Species Reactivity:	Species Independent
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Host:	Rabbit
Applications:	ELISA; the recommended starting dilution is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Serial dilutions of a BSA control, CD36 antibody, and two ADCs, 15-lipoxygenase 1 (15-LO-1) monoclonal antibody conjugated to VcMMAE (IgG VcMMAE (15-LO-1)) and CD36 monoclonal antibody conjugated to VcMMAE (IgG VcMMAE (CD36)), were plated and detected by VcMMAE Polyclonal Antibody - Biotinylated (Item No. 42295). The assay detected less than 5 ng ADC/well.

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, 04/17/2025

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

VcMMAE is a drug-linker molecule that has been used in the synthesis of antibody-drug conjugates (ADCs).^{1,2} It contains maleimide, the cathepsin B substrate valine-citrulline (Val-Cit) dipeptide, and *p*-aminobenzyl carbamate (PABC) conjugated to the antimitotic agent monomethyl auristatin E (MMAE).¹ The maleimide in VcMMAE allows it to be conjugated to cysteine residues in antibodies, which target the ADC to specific cell types. Once inside the cell, the Val-Cit dipeptide is cleaved by cathepsin B and MMAE is released.^{1,3} Formulations containing VcMMAE have been used in the treatment of diffuse large B cell lymphoma (DLBCL), Hodgkin lymphoma, systemic anaplastic large cell lymphoma, and urothelial cancer. Cayman's VcMMAE Polyclonal Antibody can be used in ELISA for the detection of ADCs containing the VcMMAE linker.

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. *Bioconjug. 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).
3. Yao, H., Jiang, F., Lu, A., *et al.* Methods to design and synthesize antibody-drug conjugates (ADCs). *Int. J. Mol. Sci.* **17**(2), 194 (2016).

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