

PRODUCT INFORMATION



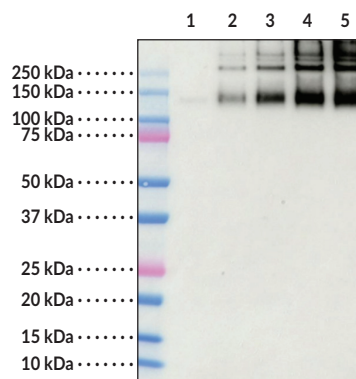
Mertansine Polyclonal Antibody - Biotinylated

Item No. 41262

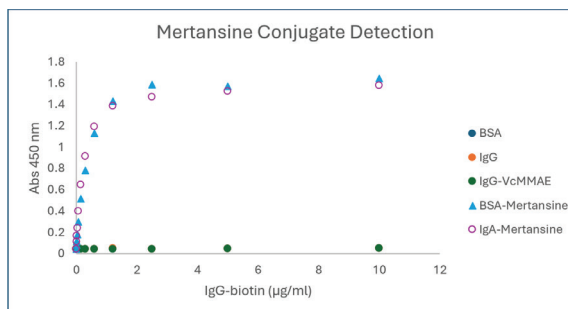
Overview and Properties

Contents:	This vial contains 100 µg of biotinylated protein A-purified rabbit IgG.
Synonyms:	DM1, Drug Maytansinoid 1, Emtansine
Immunogen:	Mertansine-KLH conjugate
Cross Reactivity:	(+) Mertansine-conjugated antibodies; (-) Albumin, VcMMAE-conjugated antibodies, Unconjugated antibodies
Species Reactivity:	Species Independent
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥2 years
Storage Buffer:	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Host:	Rabbit
Isotype:	Rabbit IgG
Applications:	ELISA and Western blot (WB); the recommended starting dilution for both applications is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

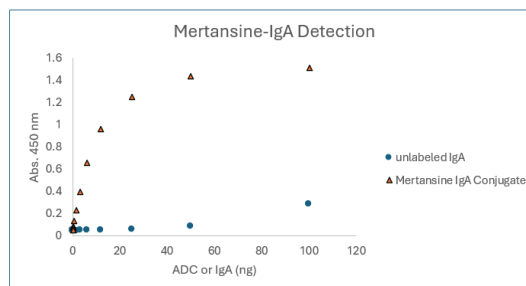
Images



Lane 1: Mouse IgA (3 µg)
Lane 2: Mouse IgA-mertansine conjugate (1 µg)
Lane 3: Mouse IgA-mertansine conjugate (2 µg)
Lane 4: Mouse IgA-mertansine conjugate (4 µg)
Lane 5: Mouse IgA-mertansine conjugate (6 µg)



Mertansine Polyclonal Antibody - Biotinylated (Item No. 41262) detected mertansine-conjugated albumin and -IgA but not unconjugated antibodies or albumin. No cross reactivity with IgG conjugated with VcMMAE (Item No. 35646).



Serial dilutions of antibody and ADC were plated and detected by Mertansine Polyclonal Antibody - Biotinylated (Item 41262). Observed a detection limit of less than Antibody - Biotinylated 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, 02/02/2026

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

Mertansine is a synthetic maytansinoid antimitotic agent.^{1,2} It disrupts mitotic spindle formation and induces cell cycle arrest and cell death in cancer cells. Mertansine has commonly been used as a payload in antibody-drug conjugates (ADCs), which have been used to selectively target and destroy cancer cells *in vitro* and *in vivo*.³⁻⁵ Formulations containing mertansine conjugated to ado-trastuzumab have been used in the treatment of HER2⁺ breast cancer. Cayman's Mertansine Polyclonal Antibody - Biotinylated can be used for ELISA and Western blot (WB) applications.

References

1. Widdison, W.C., Wilhelm, S.D., Cavanagh, E.E., *et al.* Semisynthetic maytansine analogues for the targeted treatment of cancer. *J. Med. Chem.* **49**(14), 4392-4408 (2006).
2. Lopus, M. Antibody-DM1 conjugates as cancer therapeutics. *Cancer Lett.* **307**(2), 113-115 (2011).
3. Berdeja, J.G. Lorvotuzumab mertansine: Antibody-drug-conjugate for CD56⁺ multiple myeloma. *Front. Biosci.* **19**(1), 163-170 (2014).
4. Mckertish, C.M. and Kayser, V. A novel dual-payload ADC for the treatment of HER2⁺ breast and colon cancer. *Pharmaceutics* **15**(8), 2020 (2023).
5. Huhe, M., Lou, J., Zhu, Y., *et al.* A novel antibody-drug conjugate, HcHAb18-DM1, has potent anti-tumor activity against human non-small cell lung cancer. *Biochem. Biophys. Res. Commun.* **513**(4), 1083-1091 (2019).

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