

# PRODUCT INFORMATION



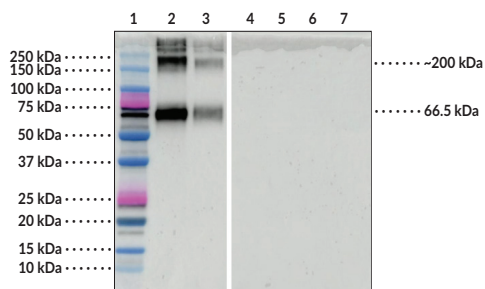
## Exatecan Monoclonal Antibody (Clone 1B2)

Item No. 44715

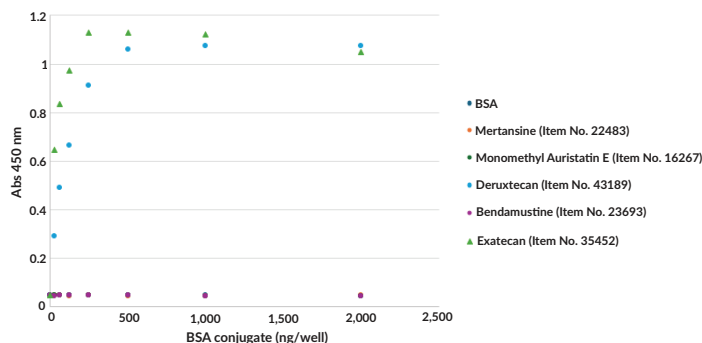
### Overview and Properties

**Contents:** This vial contains 100 µg of protein G-purified monoclonal antibody.  
**Synonym:** DX 8951f  
**Immunogen:** Exatecan-KLH  
**Cross Reactivity:** (+) Exatecan and deruxtecan; (-) Other warheads  
**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:** 1B2  
**Host:** Mouse  
**Isotype:** IgG2aK  
**Applications:** ELISA and Western blot (WB); the recommended starting dilution is 1:1,000. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Images



Lane 1: Bio-Rad Precision Plus Protein™ Dual Color Standard  
Lane 2: Exatecan (Item No. 35452)  
Lane 3: Deruxtecan (Item No. 43189)  
Lane 4: Bendamustine (Item No. 23693)  
Lane 5: Doxorubicinone (Item No. 21693)  
Lane 6: Mertansine (Item No. 22483)  
Lane 7: BSA



**ELISA.** Plates were coated with the indicated amount of BSA conjugate (or BSA alone), blocked, and probed with 1 µg./ml of the Exatecan Monoclonal Antibody (Clone 1B2) (Item No. 44715) and Goat Anti-Mouse IgG HRP (Item No. 10004302) secondary antibody.

**WB of Exatecan Monoclonal Antibody (Clone 1B2).**  
BSA conjugates (10 ng) or BSA alone were run on SDS-PAGE under reducing conditions, transferred to nitrocellulose, blocked, and probed with the Exatecan Monoclonal Antibody (Clone 1B2) (Item No. 44715) at 1 µg/ml (1:1,000) and a Goat Anti-Mouse IgG HRP (Item No. 10004302) secondary antibody.

**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, 05/27/2026

**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

---

Exatecan is a DNA topoisomerase I inhibitor that inhibits cancer cell growth *in vitro* and reduces tumor growth *in vivo*.<sup>1,2</sup> It has been used as the cytotoxic payload component of antibody-drug conjugates (ADCs), which have been used to selectively target and destroy cancer cells *in vitro* and *in vivo* in animal models.<sup>3,4</sup> Cayman's Exatecan/Deruxtecan Monoclonal Antibody (Clone 1B2) can be used for ELISA and Western blot (WB) applications.

## References

---

1. Mitsui, I., Kumazawa, E., Hirota, Y., *et al.* A new water-soluble camptothecin derivative, DX-8951f, exhibits potent antitumor activity against human tumors *in vitro* and *in vivo*. *Jpn. J. Cancer Res.* **86(8)**, 776-782 (1995).
2. Kumazawa, E. and Ochi, Y. DE-310, a novel macromolecular carrier system for the camptothecin analog DX-8951f: Potent antitumor activities in various murine tumor models. *Cancer Sci.* **95(2)**, 168-175 (2004).
3. Conilh, L., Fournet, G., Fourmaux, E., *et al.* Exatecan antibody drug conjugates based on a hydrophilic polysarcosine drug-linker platform. *Pharmaceuticals (Basel)* **14(3)**, 247 (2021).
4. Ianniello, Z., Lu, H., Quijano, E., *et al.* Harnessing AxDNA for precision exatecan delivery in cancer: A novel antibody-drug conjugate approach. *Mol. Cancer* **24(1)**, 253 (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