

PRODUCT INFORMATION



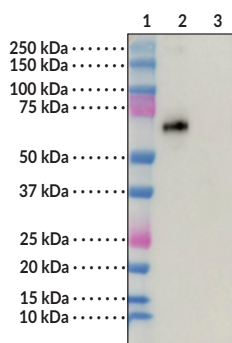
Exatecan/Dxd Polyclonal Antibody

Item No. 45399

Overview and Properties

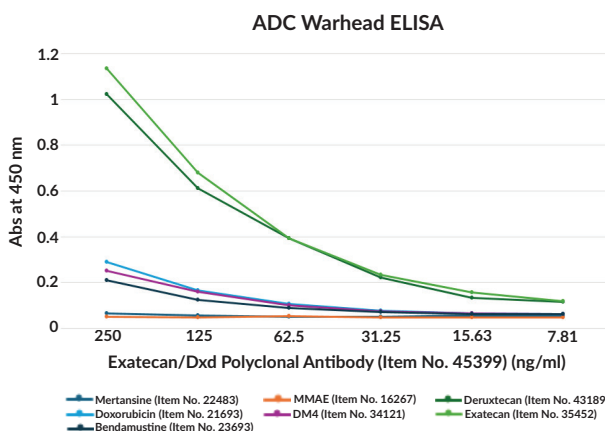
Contents: This vial contains 100 µg of protein A-purified polyclonal antibody.
Synonym: DX 8951f
Immunogen: Exatecan-KLH
Cross Reactivity: (+++) Exatecan, (++) Deruxtecan; (-) Other warheads
Species Reactivity: (+) Species independent
Form: Liquid
Storage: -20°C (as supplied)
Stability: ≥3 years
Storage Buffer: PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Host: Rabbit
Applications: ELISA and Western blot (WB); the recommended starting dilution is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Lane 1: MW Markers
Lane 2: Exatecan (Item No. 35452)
Lane 3: Bendamustine (Item No. 23693)

WB of Exatecan/Dxd Polyclonal Antibody. BSA conjugates (10 ng each) were run on an 8-16% SDS-PAGE gel under reducing conditions, then transferred to nitrocellulose and blocked with non-fat dry milk. The blot was probed with Exatecan Polyclonal Antibody (Item No. 45399) at a dilution of 1:200.



ELISA. BSA conjugates for each indicated molecule were probed with serially diluted Exatecan/Dxd Polyclonal Antibody (Item No. 45399).

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, 03/05/2026

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*.^{1,2} 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.^{3,4} Cayman's Exatecan/Dxd Polyclonal Antibody can be used for ELISA and Western blot 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