

Cyclin D1 monoclonal antibody (5D4)

Cyclins are a family of proteins which function as the regulatory subunits of cyclin/cyclin-dependent kinase (Cdk) holoenzymes, mediating entry into and progression through the cell cycle. Each cyclin is expressed in a temporal fashion during the cell cycle, partnering with a specific Cdk to activate its protein kinase activity. The most well defined mitotic cyclins are: cyclin D, which partners with cdks 4 and 6 during early G1 phase; cyclin E, which activates Cdk2 in the G1 to S phase transition; cyclin A, which utilizes Cdks 1 and 2 during progression through S phase and into G2; and cyclin B, which activates Cdk1 in M-phase promotion.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 2

[View Online »](#)

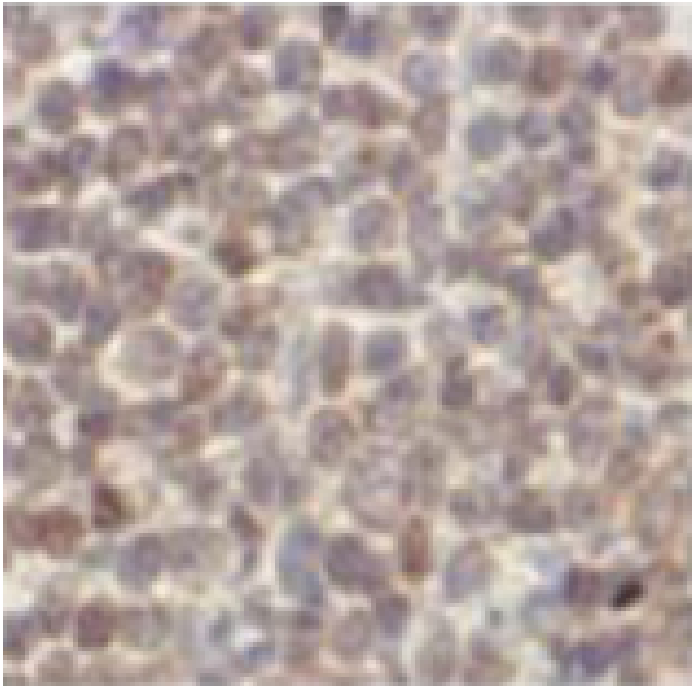
Ordering Information

[Order Online »](#)

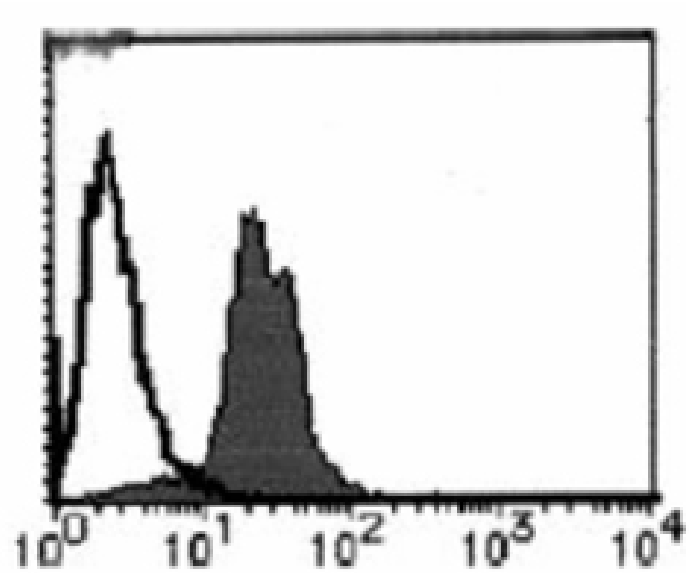
ADI-KAM-CC200-E	100µg
-----------------	-------

Manuals, SDS & CofA

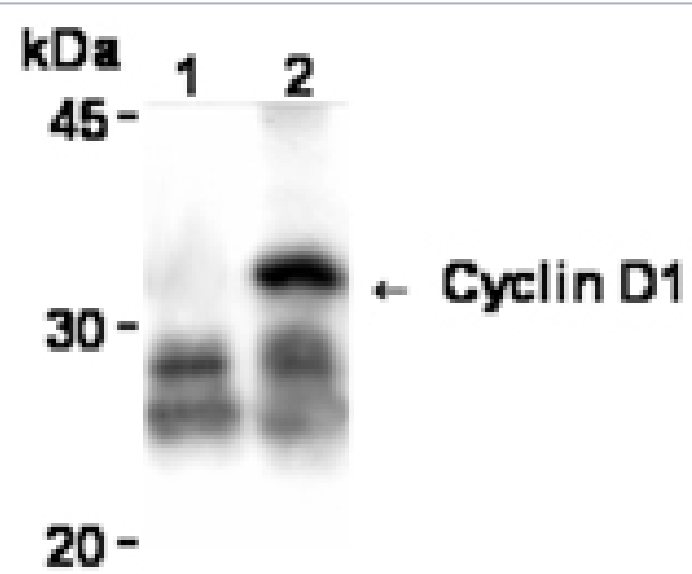
[View Online »](#)



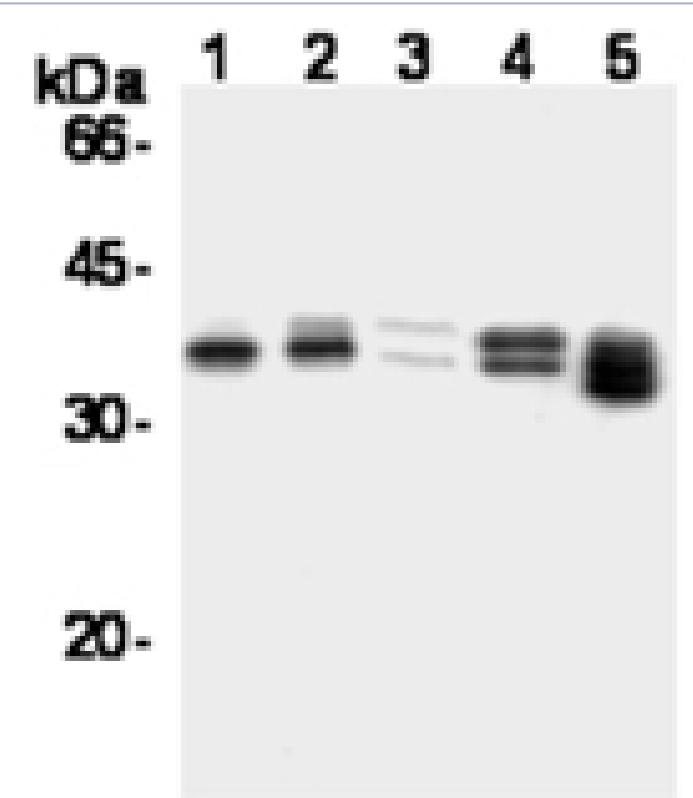
Immunohistochemistry analysis of Cyclin D1 on paraffin embedded section of human mantle cell lymphoma with Cyclin D1 mAb (5D4).



Flow cytometry analysis of Cyclin D1 expression in HeLa cells. Open histogram indicates the reaction of Isotypic control to the cells. Shaded histogram indicates the reaction of Cyclin D1 mAb (5D4) to the cells.



Immunoprecipitation analysis of cyclin D1 from ZR75-1 cells with Mouse IgG2a (1) or Cyclin D1 mAb (5D4) (2). After immunoprecipitated with the antibody, immunocomplex was resolved on SDS-PAGE and immunoBlotted with 553.



Western blot analysis of ZR- 75-1 (1), MCF7 (2), HeLa (3), WR19L (4) and NIH/3T3 (5) cells probed with Cyclin D1 mAb (5D4).

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CCND1

Application Flow Cytometry, ICC, IHC (PS), IP, WB

Application Notes Detects a band of ~36kDa by Western blot.

Clone 5D4

Formulation Liquid. In PBS, pH 7.2, containing 50% glycerol.

Host Mouse

Immunogen Recombinant human cyclin D1.

Isotype IgG2a

Purity Detail Protein A affinity purified.

Recommendation
Dilutions/Conditions Flow Cytometry (10µg/ml final concentration)Immunohistochemistry (10µg/ml, frozen/paraffin)Immunoprecipitation (1-10µg/200µg cell extract from 5×10⁶ cells)Western Blot (1µg/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human, Mouse

UniProt ID P24385

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)