

ZytoDot® 2C SPEC CCND1 Break Apart Probe

For the detection of translocations involving the CCND1 gene at 11q13.3 by chromogenic

In situ hybridization (CISH)
The ZytoDot® 2C SPEC CCND1 Break Apart Probe is designed to detect translocations involving the chromosomal region 11q13.3 harboring the CCND1 gene. Translocations involving the chromosomal region t(11;14)(q13.3;q32.3) are considered to be characteristic for mantle cell lymphomas but have also been identified in B-prolymphocytic leukemia, plasma cell myelomas, B-cell chronic lymphocytic leukemia, and in splenic lymphomas with villous lymphocytes.

Ordering Information

[Order Online »](#)

ZTV-C-3075-100	100µl
----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Cyclin D1, PRAD1

Application CISH

Quantity Sufficient for 10 tests, using 10 µl probe solution per test.

Technical Info / Product Notes For the Manual please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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