

ZytoLight® SPEC CCND1 Dual Color Break Apart Probe

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

In situ hybridization (FISH).
The ZytoLight® SPEC CCND1 Dual Color 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 other lymphoproliferative disorders, such as B-prolymphocytic leukemia, in plasma cell myelomas and B-cell chronic lymphocytic leukemia.

Ordering Information [Order Online »](#)

| | |
|----------------|-------|
| ZTV-Z-2108-50 | 50µl |
| ZTV-Z-2108-200 | 200µl |

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 FISH

Formulation Solution of salt (containing synthetic polynucleotides) in water

Quantity Sufficient for 5 tests (50 µl size) or 20 tests (200 µl size), using 10 µl probe solution per test.

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).
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