

# ZytoLight® SPEC CCND1/IGH Dual Color Dual Fusion Probe

**For the detection of the translocation t(11;14)(q13.3;q32.3) by fluorescence in situ hybridization (FISH)**

The ZytoLight® SPEC CCND1/IGH Dual Color Dual Fusion Probe is designed to detect translocation t(11;14)(q13.3;q32.3) frequently found in mantle cell lymphomas. The translocation t(11;14)(q13.3;q32.3) that involves the CCND1 and IGH gene regions is detected in up to 95% of patients with mantle cell lymphomas. However, the t(11;14) has also been identified in other lymphoproliferative disorders, such as B-prolymphocytic leukemia, and in plasma cell myelomas, B-cell chronic lymphocytic leukemia, and in splenic lymphomas with villous lymphocytes.

## Ordering Information

[Order Online »](#)

ZTV-Z-2125-50	50µl
ZTV-Z-2125-200	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, Immunoglobulin Heavy Locus, IGH@

**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](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)