

# ZytoLight® SPEC CKS1B/CDKN2C Dual Color Probe

## For the qualitative detection of amplifications/gains involving the CKS1B

The ZytoLight® SPEC CKS1B/CDKN2C Dual Color Probe is designed for the detection of gains/amplifications affecting the chromosomal region 1q21.3-q22 (CKS1B) and/or deletions of the chromosomal region 1p32.2 (CDKN2C). Chromosome abnormalities are among the most common cytogenetic findings in multiple myeloma (MM).

## Ordering Information

[Order Online »](#)

ZTV-Z-2276-50	50µl
ZTV-Z-2276-200	200µl

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Application FISH

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 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)