

# ZytoLight® SPEC CUX1/EZH2/CEN 7 Triple Color Probe

For the qualitative detection of deletions involving the human CUX1 gene in 7q22.1

and the human EZH2 gene in 7q36.1 as well as alpha satellites of chromosome 7 by fluorescence in situ hybridization (FISH)

The ZytoLight® SPEC CUX1/EZH2/CEN 7 Triple Color Probe is designed to detect losses of an entire chromosome 7 (monosomy 7) and deletions of the long arm of chromosome 7 (del(7q)). Myeloid malignancies with monosomy 7 or del(7q) respond poorly to chemotherapy and are associated with an unfavorable prognosis.

## Ordering Information

[Order Online »](#)

ZTV-Z-2214-50	50µl
---------------	------

## 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, 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)