

# ZytoLight® SPEC C19MC/19p13 Dual Color Probe

For the qualitative detection of amplifications involving the C19MC locus at 19q13.42 by

**Fluorescence in situ hybridization (FISH)**  
The ZytoLight® SPEC C19MC/19p13 Dual Color Probe is designed for the detection of amplifications of the C19MC microRNA (miRNA) cluster region located at 19q13.42. According to the 2016 WHO classification of tumors of the central nervous system, amplification of this region defines a new entity of embryonal tumors with multilayered rosettes (EMTR).

## Ordering Information

[Order Online »](#)

ZTV-Z-2274-50	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)