

# ZytoLight® SPEC NR4A3 Dual Color Break Apart Probe

For the detection of translocations involving  
the NR4A3 gene at 9q22.33-q31.1 by

Fluorescence *in situ* hybridization (FISH).

The ZytoLight® SPEC NR4A3 Dual Color Break Apart Probe is designed to detect translocations involving the chromosomal region 9q22.33-q31.1. The most frequent reciprocal translocation is t(9;22)(q22.3-q31;q12.2) found in about 70% of EMC generating a EWSR1-NR4A3 fusion gene in which the 3'-terminal part of EWSR1 is replaced by the entire NR4A3 gene. EMC is histologically characterized by a mixture of cellular and myxoid stromal components, making it difficult to distinguish it from other benign or malignant mesenchymal tumors.

## Ordering Information

[Order Online »](#)

ZTV-Z-2145-50

50µl

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Nuclear receptor subfamily 4 group A member 3, TEC, NOR1, CHN

**Application** FISH

**Quantity** Sufficient for 5 tests, 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)