

# ZytoLight® SPEC NUP214 Dual Color Break Apart Probe

For the qualitative detection of  
translocations involving the NUP214 gene at

9q34.13 by fluorescence in situ hybridization (FISH)

The ZytoLight® SPEC NUP214 Dual Color Break Apart Probe is designed for the detection of translocations involving the chromosomal region 9q34.13 harboring the NUP214 gene. Rearrangements of the NUP214 gene have been implicated in the pathogenesis of several types of hematologic malignancies, including T-cell acute lymphoblastic leukaemia (T-ALL), acute myeloid leukemia (AML), and also myelodysplastic syndrome (MDS).

## Ordering Information

[Order Online »](#)

ZTV-Z-2265-50	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)