

# ZytoLight® SPEC BCL2/IGH Dual Color Dual Fusion Probe

For the detection of the translocation  
**t(14;18)(q32.2;q21.3) by fluorescence *in situ***

**Hybridization (FISH):**  
The ZytoLight® SPEC BCL2/IGH Dual Color Dual Fusion Probe is designed to detect the translocation t(14;18)(q32.3;q21.3) affecting the BCL2 gene in the chromosomal region 18q21.33 and the IGH locus in 14q32.33. Translocations involving the BCL2 gene and the IGH gene are considered to be cytogenetic hallmarks for follicular lymphoma (FL). FL represents one of the most common non-Hodgkin lymphoma.

## Ordering Information

[Order Online »](#)

ZTV-Z-2114-50	50µl
ZTV-Z-2114-200	200µl

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** B-Cell Lymphoma 2, Immunoglobulin Heavy Locus, IGH@

**Application** FISH

**Formulation** Solution of salt (containing synthetic polynucleotides) in water

**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 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)