

# ZytoLight® SPEC IGL Dual Color Break Apart Probe

## For the qualitative detection of translocations involving the human IGL gene

The ZytoLight® SPEC IGL Dual Color Break Apart Probe is designed to detect rearrangements affecting the chromosomal region 22q11.22 at 22q11.22 by fluorescence in situ hybridization (FISH) harboring the IGL gene cluster region. The majority of translocations involve the immunoglobulin heavy chain (IGH) locus while a minor part involves the immunoglobulin light chain loci, either the kappa light chain (IGK) or the lambda light chain (IGL). IGK and IGL rearrangements resulting from the variant translocations t(2;8)(p12;q24.21) and t(8;22)(q24.21;q11.2), respectively, have been detected in up to 25% of Burkitt's Lymphoma cases.

### Ordering Information

[Order Online »](#)

ZTV-Z-2286-50	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)