

# ZytoDot® 2C SPEC BCL6 Break Apart Probe

For the detection of translocations involving  
the BCL6 gene at 3q27.3 by chromogenic *in*

**situ hybridization (CISH)**  
The ZytoDot® 2C SPEC BCL6 Break Apart Probe is designed for the detection of translocations involving the chromosomal region 3q27.3 harboring the BCL6 gene. Chromosomal rearrangements of the BCL6 gene region were found to occur in different types of non-Hodgkin lymphoma , including diffuse large B-cell lymphoma and follicular lymphoma.

## Ordering Information

[Order Online »](#)

ZTV-C-3074-100	100µl
----------------	-------

## 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 6, ZNF51, LAZ3

Application CISH

Quantity Sufficient for 10 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)