

# ZytoMation® BCL6 Dual Color Break Apart FISH Probe

## For the qualitative detection of translocations involving the human BCL6

The ZytoMation® BCL6 Dual Color Break Apart FISH Probe is designed to detect translocations involving the chromosomal region 3q27.3 harboring the BCL6 gene. Translocations involving the BCL6 gene are commonly identified in B-cell lymphomas. The most common rearrangement is the translocation t(3;14)(q27;q32.3) resulting in BCL6-IGH gene fusion. In DLBCL, the most common histologic subtype of NHL, BCL6 translocations represent one of the most frequent cytogenetic abnormality, occurring in 20% to 40% of the cases.

### Ordering Information

[Order Online »](#)

ZTV-Z-2313-5.1ML	5.1ml
------------------	-------

### 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 up to 20 tests using 240 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)