

FlexISH® BCL2/BCL6

DistinguISH Probe

For the qualitative detection of translocations involving the BCL2 gene at 18q21.33 and the BCL6 gene at 3q27.3 by fluorescence in situ hybridization (FISH)

The FlexISH® BCL2/BCL6 DistinguISH™ Probe is designed to detect rearrangements involving the chromosomal regions 18q21.33 and 3q27.3 harboring the BCL2 (B-Cell CLL/Lymphoma 2, a.k.a. PPP1R50) gene and the BCL6 (B-Cell CLL/ Lymphoma 6, a.k.a. ZNF51, LAZ3) gene, respectively. Using this probe, it is possible to simultaneously detect BCL2 and/or BCL6 rearrangements commonly found in double-hit lymphoma (DHL) or triple-hit lymphoma (THL).

Ordering Information

[Order Online »](#)

ZTV-Z-2283-50	50µl
ZTV-Z-2283-200	200µl

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application FISH

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 Manual please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com