

ZytoMation® BCL2 Dual Color Break Apart FISH Probe

For the qualitative detection of
translocations involving the human BCL2

gene at 18q21.33 by fluorescence *in situ* hybridization (FISH) on automated Bond systems

The ZytoMation® BCL2 Dual Color Break Apart FISH Probe is designed to detect translocations involving the chromosomal region 18q21.33 harboring the BCL2 gene. Translocations involving the BCL2 gene are commonly identified in B-cell lymphomas. In particular, the translocation t(14;18)(q32.3;q21.3) has been identified in about 80% of follicular lymphoma.

Ordering Information

[Order Online »](#)

ZTV-Z-2306-5.1ML	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

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