

ZytoMation® IGH Dual Color Break Apart FISH Probe

For the qualitative detection of translocations involving the human IGH

Focus at 14q32.33 by fluorescence *in situ* hybridization (FISH) on automated Bond systems

The ZytoMation® IGH Dual Color Break Apart FISH Probe is designed to detect translocations involving the chromosomal region 14q32.33 harboring the IGH locus. Rearrangements involving the IGH gene locus are considered to be cytogenetic hallmarks for non-Hodgkin lymphoma. IGH locus rearrangements have been identified in about 50% of NHLs and are associated with specific subtypes of NHLs.

Ordering Information

[Order Online »](#)

ZTV-Z-2317-5.1ML	5.1ml
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

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