

ZytoDot® 2C SPEC IGH Break Apart Probe

For the detection of translocations involving the IGH locus at 14q32.33 by chromogenic *in situ* hybridization (CISH)

The ZytoDot® 2C SPEC IGH Break Apart Probe is designed to detect translocations involving the chromosomal region 14q32.33 harboring the IGH gene. Rearrangements involving the IGH gene are considered to be cytogenetic hallmarks for non-Hodgkin lymphoma which represent 50% of all hematological malignancies.

Ordering Information

[Order Online »](#)

ZTV-C-3071-100	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 Immunoglobulin Heavy Locus, IGH@

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

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