

ZytoLight® SPEC MYC/IGH Dual Color Dual Fusion Probe

For the detection of the translocation
t(8;14)(q24;q32) by fluorescence *in situ*

Hybridization (FISH).

The ZytoLight® SPEC MYC/IGH Dual Color Dual Fusion Probe is designed to detect the translocation t(8;14)(q24;q32) affecting the MYC gene in the chromosomal region 8q24.21 and the IGH locus in 14q32.33. The most frequent translocation involving the MYC gene region t(8;14)(q24.21;q32.3) can be found in approx. 80% of the Burkitt's lymphoma cases and juxtaposes the MYC gene next to the IgH (immunoglobulin heavy chain) locus.

Ordering Information

[Order Online »](#)

ZTV-Z-2105-50	50µl
ZTV-Z-2105-200	200µl

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Protect from light.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name c-Myc, Myc Proto-oncogene, Transcription Factor p64, Immunoglobulin Heavy Locus

Application FISH

Formulation Solution of salt (containing synthetic polynucleotides) in water

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 Original Manufacturer's data sheet please [click here](#).
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