

ZytoLight® SPEC MALT1/IGH Dual Color Dual Fusion Probe

For the qualitative detection of translocations involving the human MALT1

The ZytoLight SPEC MALT1/IGH Dual Color Dual Fusion Probe (PL279) is intended to be used for the qualitative detection of the translocation t(14;18)(q32.3;q21.3) involving the human MALT1 gene at 18q21.32 and the human IGH locus at 14q32.33 in formalin-fixed, paraffin-embedded specimens by fluorescence in situ hybridization (FISH).

Ordering Information

[Order Online »](#)

ZTV-Z-2325-50	50µl
---------------	------

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

Application FISH

Quantity Sufficient for 5 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](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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