

ZytoLight® SPEC MALT1 Dual Color Break Apart Probe

For the detection of translocations involving
the MALT1 gene at 18q21.32 by fluorescence

In situ hybridization (FISH)
The ZytoLight® SPEC MALT1 Dual Color Break Apart Probe is designed to detect translocations involving the chromosomal region 18q21.32 harboring the MALT1 gene. The MALT1 gene encodes a human paracaspase and is often rearranged in MALT lymphomas accounting for 5-10% of all B-cell non-Hodgkin lymphomas.

Ordering Information

[Order Online »](#)

ZTV-Z-2196-50	50µl
ZTV-Z-2196-200	200µl

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 Mucosa associated lymphoid tissue lymphoma translocation gene 1, MLT

Application FISH

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 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