

ZytoLight® SPEC ATM/CEN 12 Dual Color Probe

For the qualitative detection of 11q22.3 deletions as well as aneusomy of

chromosome 12 by fluorescence *in situ* hybridization (FISH)

The ZytoLight® SPEC ATM/CEN 12 Dual Color Probe is designed for the detection of deletions affecting the ATM gene as well as for the enumeration of chromosome 12. In chronic lymphocytic leukemia (CLL), recurrent alterations include deletions at chromosome 13q12, 11q22.3, 17q13, and 6q21, trisomy 12 and IGH translocation. It was shown that compared to the absence of cytogenetic abnormalities, 17p13 deletion conferred the worst prognosis, followed by 11q22.3 deletion and trisomy 12.

Ordering Information

[Order Online »](#)

ZTV-Z-2296-50	50µl
---------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

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

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