

ZytoLight® SPEC TP53/ATM Dual Color Probe

For the detection of deletions affecting the human genes TP53 and ATM by fluorescence

In situ hybridization (FISH): The ZytoLight® SPEC TP53/ATM Dual Color Probe is designed for the detection of deletions affecting the genes TP53 and ATM. Chronic lymphocytic leukemia is the most common form of leukemia in Western population. TP53 gene deletions have been detected in patients with CLL, multiple myeloma, and acute myeloid leukemia. In CLL patients, allelic loss of the short arm of chromosome 17 is associated with treatment failure with alkylating agents and short survival times.

Ordering Information

[Order Online »](#)

ZTV-Z-2159-50	50µl
ZTV-Z-2159-200	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 Tumor protein 53, p53, Ataxia telangiectasia mutated

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 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