

ZytoLight® SPEC MCL1/1p12 Dual Color Probe

For the detection of the human **ML1 gene**
and **chromosome 1 specific sequences by**

Fluorescence *in situ* hybridization (FISH) The ZytoLight® SPEC MCL1/1p12 Dual Color Probe is designed for the detection of MCL1 gene amplifications. MCL1 amplifications have been reported in several human cancers including bladder, gastric, ovarian, lung, breast, melanoma, and hematologic malignancies. Overexpression of MCL1 reduces MYC-induced apoptosis in immortalized bronchial epithelial cells.

Ordering Information

[Order Online »](#)

ZTV-Z-2173-200	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 Myeloid cell leukemia 1, BCL2L3

Application FISH

Quantity Sufficient for 20 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