

ZytoLight® SPEC TP53/17q22 Dual Color Probe

For the detection of the human TP53 gene
and 17q22 specific sequences by

fluorescence *in situ* hybridization (FISH).
The ZytoLight® SPEC TP53/17q22 Dual Color Probe is designed for the detection of TP53 deletions as well as for the determination of copy number changes of the chromosomal region 17q22, harboring the MPO gene. TP53 gene deletions have been detected in patients with chronic lymphocytic leukemia, multiple myeloma, and acute myeloid leukemia.

Ordering Information

[Order Online »](#)

ZTV-Z-2198-50	50µl
---------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Tumor protein 53, p53, BCC7, LFS1, TRP53, Chromosome 17

Application FISH

Quantity Sufficient for 5 tests, 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