

ZytoLight® SPEC TP53/CEN 17 Dual Color Probe

For the detection of the human TP53 gene and alpha-satellites of chromosome 17 by

fluorescence *in situ* hybridization (FISH).

The ZytoLight® SPEC TP53/CEN 17 Dual Color Probe is designed for the detection of TP53 gene deletions observed e.g. in chronic lymphocytic leukemia (CLL). TP53 gene deletions have been detected in patients with CLL multiple myeloma (MM), acute myeloid leukemia (AML), and are also very frequent in primary solid tumors of different histological origin.

Ordering Information

[Order Online »](#)

| | |
|----------------|-------|
| ZTV-Z-2153-50 | 50µl |
| ZTV-Z-2153-200 | 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, BCC7, LFS1, TRP53, Centromere 17

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