

ZytoLight® SPEC

TERT/5q31 Dual Color

Probe

For the detection of the human TERT gene and chromosome 5 specific sequences by

fluorescence *in situ* hybridization (FISH). The ZytoLight® SPEC TERT/5q31 Dual Color Probe is designed for the detection of TERT gene amplifications and chromosomal gains found in a variety of human tumors. Chromosomal gain or amplification of the TERT gene was found in various human tumors such as lung, cervical, bladder, breast, hepatocellular and colorectal carcinomas as well as in neuroblastoma and melanoma. It was shown that TERT amplification is a poor prognostic factor in non-small cell lung cancer.

Ordering Information

[Order Online »](#)

ZTV-Z-2091-50	50µl
ZTV-Z-2091-200	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 Telomerase Reverse Transcriptase/Centromere 4

Application FISH

Formulation Solution of salt (containing synthetic polynucleotides) in water

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](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)