

ZytoLight® SPEC TERT Dual Color Break Apart Probe

For the qualitative detection of translocations involving the chromosomal

region 5p15.33 harbouring the TERT gene by fluorescence in situ hybridization (FISH)

The ZytoLight® SPEC TERT Dual Color Break Apart Probe is designed to detect translocations involving the chromosomal region 5p15.33 harboring the TERT gene. The TERT (telomerase reverse transcriptase, a.k.a. EST2, TCS1) gene encodes the reverse transcriptase component, a catalytic domain of the enzyme telomerase. Remodeling of the genomic context, due to 5p15.33 rearrangements, abrogates transcriptional silencing of the TERT gene. TERT translocations are the second common aberration found in neuroblastomas.

Ordering Information

[Order Online »](#)

ZTV-Z-2273-50	50µl
---------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application FISH

Quantity Sufficient for 5 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