

ZytoLight® SPEC TERC/CEN 3 Dual Color Probe

For the qualitative detection of amplifications involving the TERC gene at 3q26.2 by

Fluorescence in situ hybridization (FISH)
The ZytoLight® SPEC TERC/CEN 3 Dual Color Probe is designed to detect amplifications affecting the chromosomal region 3q26.2 harboring the TERC gene. 3q copy number gains including the TERC gene locus have been found in several epithelial carcinomas such as cervix carcinoma, prostate cancer, non-small cell lung cancer, and lung squamous cell carcinoma.

Ordering Information

[Order Online »](#)

ZTV-Z-2284-200	200µl
----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

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