

ZytoLight® SPEC RICTOR/5q31.1 Dual Color Probe

For the qualitative detection of amplifications involving the human RICTOR gene by

Fluorescence in situ hybridization (FISH)
The ZytoLight® SPEC RICTOR/5q31.1 Dual Color Probe is designed to detect amplifications affecting the chromosomal region 5p13.1 harboring the RICTOR. In 8-13% of non-small cell lung carcinomas (NSCLC) and small cell lung cancer (SCLC) an amplification of the RICTOR gene region has been demonstrated.

Ordering Information

[Order Online »](#)

ZTV-Z-2278-200	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