

ZytoLight® SPEC SOX2/CEN 3 Dual Color Probe

For the detection of the human SOX2 gene
and alpha-satellites of chromosome 3 by

fluorescence *in situ* hybridization (FISH).
The ZytoLight® SPEC SOX2/CEN 3 Dual Color Probe is designed for the detection of SOX2 gene amplifications frequently observed in squamous cell carcinoma of the lung, the esophagus, the oral cavity, and further organ sites. In addition, amplifications and/or overexpression were found in glioma, breast cancer, and other tumor types.

Ordering Information

[Order Online »](#)

ZTV-Z-2127-200	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 Sex Determining Region Y-Box 2, ANOP3, Centromere 3

Application FISH

Formulation Solution of salt (containing synthetic polynucleotides) in water

Quantity Sufficient for 20 tests, 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)