

ZytoLight® SPEC ROS1/CEN 6 Dual Color Probe

For the detection of the human ROS1 gene
and alpha-satellites of chromosome 6 by

fluorescence *in situ* hybridization (FISH).
The ZytoLight® SPEC ROS1/CEN 6 Dual Color Probe is designed for the detection of amplifications of the chromosomal region harboring the ROS1 gene. ROS1 has been found to undergo genetic rearrangements in a variety of human cancers including glioblastoma, cholangiocarcinoma, and non-small cell lung cancer.

Ordering Information

[Order Online »](#)

ZTV-Z-2162-200	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 c-ros Oncogene 1, MCF3, Centromere 6

Application FISH

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

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