

ZytoLight® SPEC PIK3CA/CEN 3 Dual Color Probe

For the detection of the human PIK3CA gene and alpha-satellites of chromosome 3 by

Fluorescence *in situ* hybridization (FISH).
The ZytoLight® SPEC PIK3CA/CEN 3 Dual Color Probe is designed for the detection of PIK3CA gene amplifications frequently found in a variety of human cancers. Amplifications of PIK3CA were found e.g. in cervical, ovarian, endometrial, breast, gastric, and lung cancer. In ovarian cancer as well as cervical cancer cells increased copy numbers were shown to be associated with increased expression of the gene product and PI3K activity.

Ordering Information

[Order Online »](#)

ZTV-Z-2140-200	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 Phosphatidylinositol-4,5-Bisphosphate 3-Kinase, Catalytic Subunit Alpha, PI3K-alpha, 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)