

ZytoLight® Aneusomy Probe Set

For the qualitative detection of alpha-satellites of chromosome X and the classical satellites III of chromosome Y as well as alpha satellites of chromosome 18 and

Human chromosome 13 and 21 specific sequences by fluorescence in situ hybridization (FISH)

The ZytoLight® Aneuploidy Panel XY and 13/18/21 is designed for enumeration of the chromosomes 13, 18, 21, X, and Y. Trisomies of the autosomes 13, 18, or 21 (Down Syndrome) are common genomic aberrations. Aberrant numbers of the gonosomes X and Y are resulting in disorders of sex development.

Ordering Information

[Order Online »](#)

ZTV-Z-2104-5	5 tests
ZTV-Z-2104-20	20 tests

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Protect from light.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application FISH

Contents ZytoLight CEN X/Y Dual Color Probe (Prod. No. [ZTV-Z-2016](#))
ZytoLight SPEC 13/CEN 18/SPEC 21 Triple Color Probe (Prod. No. [ZTV-Z-2095](#))

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)