

ZytoLight® CEN 10 Probe (Alpha Sat. D10Z1)

For the detection of human alpha-satellites of chromosome 12 by fluorescence *in situ* hybridization (FISH).

Ordering Information

[Order Online »](#)

ZTV-Z-2079-200	200µl
----------------	-------

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

Alternative Name Centromere 10

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

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