

ZytoLight® SPEC

D13S319/13q34/CEN

12 Triple Color Probe

For the detection of the D13S319 locus,
human 13q34 specific sequences, and alpha-

satellites of chromosome 12 by fluorescence
in situ hybridization (FISH).

The ZytoLight® SPEC D13S319/13q34/ CEN 12 Triple Color Probe is designed for the detection of D13S319 deletions as well as for the enumeration of chromosome 12. Chronic lymphocytic leukemia is the most common form of leukemia in Western population. Trisomy 12 represents another frequent chromosomal aberration in CLL, detected in about 20% of CLL cases. Trisomy 12 as single aberration is associated with an intermediate prognostic outcome.

Ordering Information

[Order Online »](#)

ZTV-Z-2160-50	50µl
ZTV-Z-2160-200	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 Tumor protein 53, p53, Ataxia telangiectasia mutated, Centromere 12

Application FISH

Quantity Sufficient for 5 tests (50 µl size) or 20 tests (200 µl size), 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