

ZytoLight® SPEC

D13S319/13q34 Dual Color Probe

For the qualitative detection of deletions involving the human D13S319 region and human 13q34 specific sequences by fluorescence in situ hybridization (FISH)

The ZytoLight® SPEC D13S319/13q34 Dual Color Probe is designed for the detection of deletions on the long arm of chromosome 13. Chronic lymphocytic leukemia (CLL) is the most common form of leukemia in the Western population. The most frequent aberration in CLL is the deletion of 13q14.2 which involves the D13S319 locus and which is associated with a favorable prognosis if occurring as the sole genetic aberration.

Ordering Information

[Order Online »](#)

ZTV-Z-2280-50-R	50µl
ZTV-Z-2280-200-R	200µl

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

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 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