

ZytoLight® SPEC MYB/CEN 6 Dual Color Probe

For the qualitative detection of deletions involving the human MYB gene and the

detection of alpha satellites of chromosome 6 by fluorescence in situ hybridization (FISH)
The ZytoLight® SPEC MYB/CEN 6 Dual Color Probe is designed to detect deletions affecting the chromosomal region 6q23.3 harboring the MYB gene. 3-10% of chronic lymphocytic leukemia (CLL) cases have been shown to harbor structural aberrations in the chromosomal region of 6q, one of them is 6q23.

Ordering Information

[Order Online »](#)

| | |
|---------------|------|
| ZTV-Z-2281-50 | 50µl |
|---------------|------|

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, 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