

ZytoLight® SPEC MYB Dual Color Break Apart Probe

For the detection of translocations involving
the MYB gene at 6q23.3 by fluorescence *in*

situ hybridization (FISH).
The ZytoLight® SPEC MYB Dual Color Break Apart Probe is designed to detect translocations involving the chromosomal region 6q23.3 harboring the MYB gene. Rearrangements of the MYC locus have been detected in adenoid cystic carcinoma as well as T-ALL.

Ordering Information

[Order Online »](#)

ZTV-Z-2143-50	50µl
ZTV-Z-2143-200	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 v-myb avian myeloblastosis viral oncogene homolog, c-Myb

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