

ZytoLight® SPEC GATA2/MECOM Dual Color Dual Fusion Probe

The ZytoLight® SPEC GATA2/MECOM Dual Color Dual Fusion Probe is designed to detect the inversion $inv(3)(q21;q26.2)$ and the translocation $t(3;3)(q21;q26.2)$ both affecting the GATA2 gene in the chromosomal region 3q26.2. $inv(3)$, $t(3;3)$, and less commonly $ins(3;3)$ occur in acute myeloid leukemia (AML) and are also observed in MDS and dysplastic syndromes and in the blastic phase of chronic myeloid leucemia (CML).

Ordering Information

[Order Online »](#)

ZTV-Z-2287-50	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