

ZytoLight® SPEC FGFR3/IGH Dual Color Dual Fusion Probe

For the qualitative detection of the
translocation **t(4;14)(p16;q32.3)** by

Fluorescence in situ hybridization (FISH)
The ZytoLight® SPEC FGFR3/IGH Dual Color Dual Fusion Probe is designed to detect the translocation t(4;14)(16.3;q32.3) affecting the FGFR3 gene in the chromosomal region 4p16.3 and IGH locus in 14q32.33. FGFR3/IGH translocations are observed in approximately 15-20% of patients with multiple myeloma (MM).

Ordering Information

[Order Online »](#)

ZTV-Z-2282-50	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