

FlexISH® MYC/IGH TriCheck™ Probe

For the qualitative detection of human MYC rearrangements with and without participation of the human IGH locus by fluorescence *in situ* hybridization (FISH)

The FlexISH® MYC/IGH TriCheck™ Probe is designed to detect the translocation t(8;14)(q24.21;q32.3) affecting the MYC gene in the chromosomal region 8q24.21 and the IGH locus in 14q32.33. Moreover, using this probe it is possible to discriminate between MYC-IGH translocations and MYC translocations involving fusion partners other than IGH. The MYC proto-oncogene encodes a transcription factor essential for cell growth and proliferation and is broadly implicated in tumorigenesis.

Ordering Information

[Order Online »](#)

ZTV-Z-2293-50	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