

FlexISH® RET/KIF5B TriCheck Probe

For the qualitative detection of human RET-KIF5B rearrangements by fluorescence in situ hybridization (FISH)

The FlexISH® RET/KIF5B TriCheck™ Probe is designed to detect inversions involving the chromosomal regions 10q11.21 and 10q11.22 harboring the RET and KIF5B genes, respectively. Moreover, using this probe it is possible to discriminate between KIF5B-RET inversions and RET translocations involving fusion partners other than KIF5B. RET rearrangement, including inversions and translocations, are found in NSCLC with an incidence of 1-2%.

Ordering Information

[Order Online »](#)

ZTV-Z-2269-200	200µl
----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

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 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