

FlexISH®

NTRK1/NTRK3

DistinguISH™ Probe

For the detection of translocations involving the human NTRK1 gene at 1q23.1 and the human NTRK3 gene at 15q25.3 by fluorescence in situ hybridization (FISH)

The FlexISH® NTRK1/NTRK3 DistinguISH™ Probe is designed to detect rearrangements affecting the chromosomal region 1q23.1 and 15q25.3 harboring the NTRK1 and NTRK3 gene region, respectively. The treatment of patients with NTRK fusion-positive cancers with a NTRK inhibitor is associated with high response rates, regardless of NTRK gene, fusion partner, and tumor type.

Ordering Information

[Order Online »](#)

ZTV-Z-2314-50	50µl
ZTV-Z-2314-200	200µl

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

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