

ZytoLight® SPEC NTRK2 Dual Color Break Apart Probe

For the detection of translocations involving the NTRK2 gene at 9q21.33 by fluorescence

In situ hybridization (FISH)
The ZytoLight® SPEC NTRK2 Dual Color Break Apart Probe is designed to detect translocations involving the chromosomal region 9q21.33 harboring the NTRK2 gene. NTRK2 is a tyrosine kinase and plays a key role in central and peripheral nervous system development as well as in cell survival. Translocations affecting the NTRK2 gene have been reported in several cancer types, including glioblastomas, pilocytic astrocytomas, head and neck squamous cell carcinoma, and lung adenocarcinoma. The treatment of patients with NTRK1, 2, or 3 fusion-positive cancers with an NTRK inhibitor, such as the FDA-approved drugs larotrectinib or entrectinib, is associated with high response rates, regardless of NTRK gene, fusion partner, and tumor type.

Ordering Information

[Order Online »](#)

ZTV-Z-2205-50	50µl
ZTV-Z-2205-200	200µl

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Neurotrophic receptor tyrosine kinase 2, TRKB

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