

ZytoLight® SPEC KIF5B Dual Color Break Apart Probe

For the detection of translocations involving
the KIF5B gene at 10p11.22 by fluorescence

In situ hybridization (ISH):
The ZytoLight® SPEC KIF5B Dual Color Break Apart Probe is designed to detect translocations involving the chromosomal region 10p11.22 harboring the KIF5B gene. About 5% of all non-small cell lung cancer cases are positive for the ALK-EML4 fusion as a result of an inversion in chromosome 2.

Ordering Information

[Order Online »](#)

ZTV-Z-2131-50	50µl
---------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Kinesin Family Member 5B

Application FISH

Formulation Solution of salt (containing synthetic polynucleotides) in water

Quantity Sufficient for 5 tests, using 10 µl probe solution per test.

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).
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