

ZytoLight® SPEC

ALK/2q11 Dual Color

Probe

For the detection of the human ALK gene and chromosome 2 specific sequences by

Fluorescence *in situ* hybridization (FISH).

The ZytoLight® SPEC ALK/2q11 Dual Color Probe is designed for detection of amplifications of the chromosomal region harboring the ALK gene. In addition to ALK gene fusions, ALK has also shown to be amplified and show copy number gains in a variety of tumors including non-small cell lung cancer and alveolar rhabdomyosarcoma. In colorectal cancer, ALK amplification was correlated with nodal status suggesting that ALK amplified tumors are also a frequent genetic event in the tumorigenesis of neuroblastomas and were found to result in high ALK expression correlating with an unfavorable neuroblastoma phenotype.

Ordering Information

[Order Online »](#)

ZTV-Z-2161-200	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 Anaplastic Lymphoma Receptor Tyrosine Kinase, CD246, Chromosome 2

Application FISH

Quantity Sufficient for 20 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