

FlexISH® ALK/ROS1 DistinguISH™ Probe

For the detection of translocations involving the ranslocations involving the ALK gene at 2p23.1-p23.2 and the ROS1 gene at 6q22.12 by fluorescence *in situ* hybridization (FISH)

The FlexISH® ALK/ROS1 DistinguISH™ Probe is designed to detect rearrangements involving the ALK (anaplastic lymphoma receptor tyrosine kinase) and ROS1 (c-ros oncogene 1) gene, respectively. Using this probe, it is possible to simultaneously detect ALK and ROS1 rearrangements and, additionally, to discriminate between possible aberrations affecting these chromosomal regions.

Ordering Information

[Order Online »](#)

ZTV-Z-2203-50	50µl
ZTV-Z-2203-200	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, c-ros oncogene 1

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