

# ZytoMation® RET Dual Color Break Apart FISH Probe

For the qualitative detection of  
translocations involving the human RET

gene at 10q11.21 by fluorescence *in situ* hybridization (FISH) on automated Bond systems

The ZytoMation® RET Dual Color Break Apart FISH Probe is designed to detect translocations involving the chromosomal region 10q11.21 harboring the RET gene. RET encodes a tyrosine kinase receptor. RET translocations are responsible for 1-2% of non-squamous NSCLCs. Similarly to ALK and ROS1, they are more characteristic for young non-smokers and females.

## Ordering Information

[Order Online »](#)

ZTV-Z-2316-5.1ML	5.1ml
------------------	-------

## 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 up to 20 tests using 240 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](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)