

# ZytoDot® 2C SPEC RET Break Apart Probe

For the detection of translocations involving the RET gene at 10q11.21 by chromogenic *in situ* hybridization (CISH)

The ZytoDot® 2C SPEC RET Break Apart Probe is designed to detect translocations involving the chromosomal region 10q11.21 harboring the RET gene. Translocations involving RET were first described in papillary thyroid carcinoma where somatic rearrangements result in the fusion of its TK catalytic domain with an N-terminal dimerization domain encoded by various fusion partner genes.

## Ordering Information

[Order Online »](#)

ZTV-C-3064-100	100µl
----------------	-------

## 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** Rearranged during transfection proto-oncogene

**Application** CISH

**Quantity** Sufficient for 10 tests (100 µl size) or 40 tests (400 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](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)