

# ZytoLight® SPEC RREB1/MYB/CEN 6 Triple Color Probe

For the detection of the human RREB1 gene,  
the human MYB gene, and alpha-satellites of

chromosome 6 by fluorescence *in situ* hybridization (FISH)

The ZytoLight® SPEC RREB1/MYB/CEN 6 Triple Color Probe is designed for the detection of copy number changes of the chromosomal regions harboring the RREB1 and the MYB gene, respectively. Overexpression of the RREB1 protein was detected in prostate cancer and in a medullary thyroid cancer cell line. MYB has been found to be amplified in a variety of human cancers. In pancreatic cancer, MYB amplification was mainly found in advanced and metastatic tumors.

## Ordering Information

[Order Online »](#)

ZTV-Z-2152-50	50µl
ZTV-Z-2152-200	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** Ras responsive element binding protein 1, HNT, v-myb avian myeloblastosis viral oncogene homolog, c-myb, Centromere 6

**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](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)