

# ZytoLight® SPEC ERBB2/CEN 17 Dual Color Probe

For the detection of human ERBB2 gene and  
alpha-satellites of chromosome 17 by

**Fluorescence *in situ* hybridization (FISH).**  
The ZytoLight® SPEC ERBB2/CEN 17 Dual Color Probe is designed for  
the detection of ERBB2 (a.k.a HER2) gene amplification frequently  
observed in solid malignant neoplasms e.g. breast cancer samples.

## Ordering Information

[Order Online »](#)

ZTV-Z-2015-50	50µl
ZTV-Z-2015-200	200µl

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Handling** Protect from light.

**Long Term Storage** +4°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** c-erbB-2, HER-2, Centromere 17

**Application** FISH

**Formulation** Solution of salt (containing synthetic polynucleotides) in water

**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 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](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)