

# ZytoLight® SPEC FGFR2/CEN 10 Dual Color Probe

For the detection of the human FGFR2 gene  
and alpha-satellites of chromosome 10 by

**Fluorescence *in situ* Hybridization (FISH).**  
The ZytoLight® SPEC FGFR2/CEN 10 Dual Color Probe is designed for the detection of FGFR2 gene amplifications frequently observed in breast cancer as well as in gastric cancer. Moreover, high FGFR2 expression is correlated with poor overall survival and poor disease-free survival rates in breast cancer patients.

## Ordering Information

[Order Online »](#)

ZTV-Z-2122-200	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** Fibroblast Growth Factor Receptor 1/Centromere 10

**Application** FISH

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

**Quantity** Sufficient for 20 tests, 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)