

# ZytoLight® SPEC VEGFA/CEN 6 Dual Color Probe

For the detection of the human VEGFA gene  
and alpha-satellites of chromosome 6 by

**Fluorescence *in situ* hybridization (FISH)**

The ZytoLight® SPEC VEGFA/CEN 6 Dual Color Probe is designed for the detection of amplifications involving the chromosomal region 6p21.1 harboring the VEGFA gene. Amplification of the VEGFA gene locus was found in several types of malignancy, such as osteosarcoma, hepatocellular carcinoma, and colorectal cancers.

## Ordering Information

[Order Online »](#)

ZTV-Z-2195-200	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 Vascular endothelial growth factor A, Centromere 6

Application FISH

Quantity Sufficient for 20 tests, 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)