

# ZytoLight® SPEC COL1A1/PDGFB Dual Color Dual Fusion Probe

**For the detection of the translocation t(17;22)(q21.3;q13.1) by fluorescence in situ hybridization (FISH)**

The ZytoLight® SPEC COL1A1/PDGFB Dual Color Dual Fusion Probe is designed for the detection of the specific translocations involving the chromosomal region 17q21.33 harboring the COL1A1 gene, and the chromosomal region 22q13.1, harboring the PDGFB gene. The reciprocal translocations involving t(17;22)(q21.3;q13.1) are characteristic for dermatofibrosarcoma protuberans (DFSP) patients. DFSP is a highly recurrent, infiltrative skin tumor of intermediate malignancy.

## Ordering Information

[Order Online »](#)

ZTV-Z-2116-50	50µl
ZTV-Z-2116-200	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** Collagen Type I Alpha 1/Platelet-Derived Growth Factor Beta Polypeptide

**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)