

# ZytoLight® SPEC FUS Dual Color Break Apart Probe

For the detection of translocations involving  
the FUS gene at 16p11.2 by fluorescence *in*

situ hybridization (FISH).

The ZytoLight® SPEC FUS Dual Color Break Apart Probe is designed to detect translocations involving the chromosomal region 16p11.2 harboring the FUS gene. FUS gene rearrangements have been shown to be involved in both solid tumors and leukemias. Occurring in over 90% of myxoid liposarcomas, the FUS-DDIT3 fusion protein is regarded as being consequential for the development of myxoid liposarcomas.

## Ordering Information

[Order Online »](#)

|               |      |
|---------------|------|
| ZTV-Z-2130-50 | 50µl |
|---------------|------|

## 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** Fused In Sarcoma, ER81 TLS, HNRNPP2

**Application** FISH

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

**Quantity** Sufficient for 5 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)