

ZytoLight® SPEC PHF1 Dual Color Break Apart Probe

For the qualitative detection of translocations involving the human PHF1

gene at 6p21.32 by fluorescence in situ hybridization (FISH)

The ZytoLight® SPEC PHF1 Dual Color Break Apart Probe is designed for the detection of translocations involving the chromosomal region 6p21.32 harboring the PHF1 gene. Endometrial stromal tumors (ESTs) are the second most common pure mesenchymal tumors of the uterus. ESTs may pose diagnostic challenges particularly when they exhibit variant histologic appearances.

Ordering Information

[Order Online »](#)

ZTV-Z-2215-50	50µl
---------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application FISH

Quantity Sufficient for 5 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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com