

ZytoLight® SPEC JAZF1 Dual Color Break Apart Probe

For the detection of translocations involving
the JAZF1 gene at 7p15.1-p15.2 by

Fluorescence *in situ* hybridization (FISH).

The ZytoLight® SPEC JAZF1 Dual Color Break Apart Probe is designed to detect translocations involving the chromosomal region 7p15.1-p15.2 harboring the JAZF1 gene. In 25-30% of endometrial stromal sarcoma the JAZF1 gene is disrupted by another translocation t(6;7) where the first zinc finger domain of JAZF1 is fused to both zinc finger domains of the PHF1 gene at 6p21.32.

Ordering Information

[Order Online »](#)

| | |
|---------------|------|
| ZTV-Z-2132-50 | 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 Juxtaposed With Another Zinc Finger 1

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

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