

ZytoLight SPEC PLAG1 Dual Color Break Apart Probe

For the qualitative detection of rearrangements involving the human PLAG1 gene at 8q12.1 by fluorescence in situ hybridization (FISH)

The ZytoLight SPEC PLAG1 Dual Color Break Apart Probe (PL280) is intended to be used for the qualitative detection of rearrangements involving the human PLAG1 gene at 8q12.1 in formalin-fixed, paraffin-embedded specimens by fluorescence in situ hybridization (FISH). The probe is intended to be used in combination with the ZytoLight FISH-Tissue Implementation Kit (Prod. No. Z-2028-5/-20).

Ordering Information

[Order Online »](#)

ZTV-Z-2326-50	50µl
---------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Protect from light.

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