

ZytoLight® SPEC RB1/13q34 Dual Color Probe

For the qualitative detection of human RB1 gene deletions and human 13q34 specific

sequences by fluorescence in situ hybridization (FISH)

The ZytoLight SPEC RB1/13q34 Dual Color Probe (PL278) is intended to be used for the qualitative detection of human RB1 gene deletions as well as the detection of human 13q34 specific sequences in cytologic or formalin-fixed, paraffin-embedded specimens by fluorescence in situ hybridization (FISH). The probe is intended to be used in combination with ZytoLight FISH Implementation Kits (Prod. No. Z-2028-5/-20, or Z-2099-20).

Ordering Information

[Order Online »](#)

ZTV-Z-2324-50	50µl
ZTV-Z-2324-200	200µl

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

Alternative Name RB transcriptional corepressor 1

Application FISH

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 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