

ZytoLight® SPEC NUP98 Dual Color Break Apart Probe

For the qualitative detection of translocations involving the human NUP98

gene at 11p15.4 by fluorescence in situ hybridization (FISH)

The ZytoLight® SPEC NUP98 Dual Color Break Apart Probe is designed for the detection of translocations involving the chromosomal region 11p15.4 harboring the NUP98 gene. Rearrangements of the NUP98 gene have been implicated in the pathogenesis of several types of hematologic malignancies, including de novo and therapy-related acute myeloid leukaemia (AML), and also myelodysplastic syndrome (MDS), chronic myelogenous leukaemia (CML), and T-cell acute lymphoblastic leukaemia (T-ALL).

Ordering Information

[Order Online »](#)

ZTV-Z-2266-50	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