

ZytoLight® SPEC NUTM1 Dual Color Break Apart Probe

For the detection of translocations involving
the NUTM1 gene at 15q14 by fluorescence *in*

situ hybridization (FISH)
The ZytoLight® SPEC NUTM1 Dual Color Break Apart Probe is designed to detect translocations involving the chromosomal region 15q14 harboring the NUTM1 gene. NUT midline carcinoma (NMC) is a rare and aggressive form of squamous cell carcinoma that arises mainly in the head, neck, or mediastinum. NMCs may be indistinguishable from more common squamous cell carcinomas and are thus an underdiagnosed entity. Therefore, the diagnosis of NMC depends on the confirmation of NUTM1 rearrangement.

Ordering Information

[Order Online »](#)

ZTV-Z-2208-200	200µl
----------------	-------

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 NUT midline carcinoma family member 1, NUT

Application FISH

Quantity Sufficient for 20 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