

ZytoLight® SPEC NRG1/CD74 TriCheck Probe

For the detection of NRG1-CD74
rearrangements by fluorescence *in situ*
hybridization (FISH)

The ZytoLight® SPEC NRG1/CD74 TriCheck™ Probe is designed to detect translocations involving the NRG1 gene and the CD74 gene. Using this probe it is possible to discriminate between CD74-NRG1 fusions and translocations affecting NRG1, but not CD74, such as SLC3A2-NRG1 or VAMP2-NRG1 fusions detected in various tumors.

Ordering Information

[Order Online »](#)

ZTV-Z-2194-200	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 Neuregulin 1, HKL, GGF

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