

ZytoLight® SPEC ERBB2/TOP2A/CEN 17 Triple Color Probe

For the detection of the human ERBB2 gene,
the human TOP2A gene and alpha-satellites

of chromosome 17 by fluorescence *in situ*
hybridization (FISH).

The ZytoLight® SPEC ERBB2/TOP2A/CEN 17 Triple Color Probe is designed for the simultaneous detection of ERBB2 and TOP2A gene status. The TOP2A gene is frequently either co-amplified or deleted in ERBB2 positive breast cancer cases. TOP2A functions as the target for several anticancer agents, e.g. anthracyclines.

Ordering Information

[Order Online »](#)

ZTV-Z-2093-50	50µl
ZTV-Z-2093-200	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 c-erbB2, HER-2, Topoisomerase (DNA) II alpha, Centromere 17

Application FISH

Formulation Solution of salt (containing synthetic polynucleotides) in water

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 Original Manufacturer's data sheet please [click here](#).
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