

ZytoLight® SPEC ERBB2/CEN 17 Dual Color Probe Kit

For the detection of human ERBB2 gene and alpha-satellites of chromosome 17 by

Fluorescence *in situ* hybridization (FISH). The ZytoLight® SPEC ERBB2/CEN 17 Dual Color Probe is designed for the detection of ERBB2 (a.k.a HER2) gene amplification frequently observed in solid malignant neoplasms e.g. breast cancer samples.

Ordering Information

[Order Online »](#)

ZTV-Z-2020-5	5 tests
ZTV-Z-2020-20	20 tests

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 c-erbB-2, HER-2, Centromere 17

Application FISH

Contents Heat Pretreatment Solution Citric; Pepsin Solution; Wash Buffer SSC; Probe; 25x Wash Buffer A; DAPI/Antifade-Solution, good for 5 or 20 tests.

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](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)