

ZytoDot® 2C SPEC ERBB2/CEN 17 Probe (Digoxigenin/DNP)

For the qualitative detection of human ERBB2 gene amplifications and alpha

**satellites of chromosome 17 by chromogenic
in situ hybridization (CISH)**
For the detection of the human ERBB2 gene and alpha-satellites of
chromosome 17 by chromogenic *in situ* hybridization (CISH).

The ZytoDot 2C SPEC ERBB2/CEN 17 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. Amplification of the proto-oncogene ERBB2, observed in approximately 20% of all breast cancer samples, has been correlated with a poor prognosis of the disease.

In addition, detection of ERBB2 amplification by CISH may help in selecting patients eligible for ERBB2-targeted therapy.

Ordering Information

[Order Online »](#)

ZTV-C-3032-100	100µl
ZTV-C-3032-400	400µl

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, NEU, Centromere 17

Application CISH

Formulation Solution of salt (containing synthetic polynucleotides) in water

Quantity Sufficient for 10 tests (100 µl size) or 40 tests (400 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