

ZytoDot® SPEC

ERBB2 Probe

(Digoxigenin)

For the detection of the human ERBB2 gene by chromogenic *in situ* hybridization (CISH).

The ZytoDot® SPEC ERBB2 Probe is designed for the detection of ERBB2 gene amplification, frequently observed in solid malignant neoplasms, in tissue sections or cell 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. Similar results have been obtained for a variety of other malignant neoplasms e.g. ovarian cancer, stomach cancer, and carcinomas of the salivary gland.

Ordering Information

[Order Online »](#)

ZTV-C-3001-400	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, HER2, NEU

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