

ZytoDot® SPEC CMYC Probe (Digoxigenin)

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

The ZytoDot® SPEC MYC Probe is designed for the detection of MYC gene amplification frequently observed in malignant tumors e.g. breast and endometrial cancer. Amplification of the chromosomal MYC gene region has been detected in many types of malignant neoplasms e.g. breast, lung, head, colon, kidney, neck, ovary, bladder, and endometrial cancers.

Ordering Information

[Order Online »](#)

ZTV-C-3013-400	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-Myc, Myc Proto-oncogene, Transcription Factor p64

Application CISH

Formulation Solution of salt (containing synthetic polynucleotides) in water

Quantity Sufficient for 40 tests, 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