

ZytoDot® 2C SPEC EGFR/CEN 7 Probe (Digoxigenin/DNP)

For the detection of the human EGFR gene
and alpha-satellites of chromosome 7 by

chromogenic *in situ* hybridization (CISH). The ZytoDot® 2C SPEC EGFR/CEN 7 Probe is designed for the simultaneous detection of EGFR and centromere 7 in tissue sections and cell samples. Overexpression of EGFR has been shown in a number of tumor entities and is associated with poor prognosis.

Ordering Information

[Order Online »](#)

ZTV-C-3033-100	100µl
ZTV-C-3033-400	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 EGF Receptor, Epidermal Growth Factor Receptor, HER-1, Centromere 7

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