

ZytoDot® SPEC ERBB2 Probe Kit (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 (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.

Ordering Information

[Order Online »](#)

ZTV-C-3003-40	40 tests
---------------	----------

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

Contents Heat Pretreatment Solution EDTA; Pepsin Solution; Probe; Wash Buffer SSC; PBS/Tween; Blocking Solution; Mouse-anti-DIG; Anti-Mouse-HRP-Polymer; DAB Solution A; DAB Solution B; Mayer's Hematoxylin Solution; Mounting Solution (alcoholic); Control Slides, good for 10 or 40 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

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