

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

For the qualitative detection of human  
ERBB2 gene amplifications 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.

## Ordering Information

[Order Online »](#)

|               |          |
|---------------|----------|
| ZTV-C-3022-10 | 10 tests |
| ZTV-C-3022-40 | 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, Centromere 17

**Application** CISH

**Contents** Heat Pretreatment Solution EDTA; Pepsin Solution; Probe; Wash Buffer SSC; 20x Wash Buffer TBS; Anti-Dig/DNP-Mix; HRP/AP-Polymer-Mix; AP-Red Solution A; AP-Red Solution B; HRP-Green Solution A; HRP-Green Solution B; Nuclear Blue 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](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)