

# ZytoDot® 2C CISH Implementation Kit (Digoxigenin/DNP)

For chromogenic *in situ* hybridization (CISH)  
using any *ZytoDot* CISH Probe

## Ordering Information

[Order Online »](#)

ZTV-C-3044-10	10 tests
ZTV-C-3044-40	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

**Contents** Heat Pretreatment Solution EDTA; Pepsin Solution; 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), good for 10 or 40 Tests.

**Technical Info / Product Notes** 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)