

# Anti-DIG/Anti-DNP Mix

For use in chromogenic *in situ* hybridization (CISH) procedures

Antibodies for the detection of digoxigenin (DIG)-labeled and dinitrophenyl (DNP)-labeled probes in either formalin-fixed, paraffin-embedded tissue or cell samples by chromogenic *in situ* hybridization (CISH). Component of ZytoDot 2C CISH Polymer Detection Kit (Prod. No. [ZTV-C-3028](#)).

## Ordering Information

[Order Online »](#)

ZTV-AB-0014-4	4ml
---------------	-----

## 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** Antibodies to Digoxigenin and Dinitrophenyl

**Formulation** Solution of salt (containing synthetic polypeptides) in water

**Immunogen** Digoxigenin/DNP

**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)