

ZytoFast® PLUS CISH Implementation Kit AP-Permanent Red

For chromogenic *in situ* hybridization (CISH)
using any digoxigenin-labeled *ZytoFast* CISH
Probe

Ordering Information

[Order Online »](#)

ZTV-T-1151-40	40 tests
---------------	----------

Manuals, SDS & CofA

[View Online »](#)

CISH Accessories

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Contents DNA (+)-Control Probe, 0.1 ml; DNA (-)-Control Probe, 0.1 ml; RNA (+)-Control Probe, 0.1 ml; RNA (-)-Control Probe, 0.1 ml; Heat Pretreatment Solution EDTA, 500 ml; Pepsin Solution, 4 ml; 20x Wash Buffer TBS, 4×50 ml; Rabbit-anti-DIG, 4 ml; Anti-Rabbit-AP-Polymer, 4 ml; Permanent Red Solution A, 0.2 ml; Permanent Red Solution B, 14 ml; Mayer's Hematoxylin Solution, 20 ml; Mounting Solution (alcoholic), 4 ml

Technical Info / Product Notes 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