

Anti-Biotin/DIG-Mix

For use in chromogenic in situ hybridization (CISH) procedures

The Anti-Biotin/DIG-Mix (AB15) is intended to be used for detection steps in chromogenic in situ hybridization (CISH) applications on formalin-fixed, paraffin-embedded specimens. The Anti-Biotin/DIG-Mix is intended to be used in combination with either the ZytoFast human Ig-kappa/Ig-lambda CISH Kit (Prod. No. T-1005-40), or the ZytoFast human Ig-kappa/Ig-lambda Permanent CISH Kit (Prod. No. T-1105-40).

Ordering Information

[Order Online »](#)

ZTV-AB-0015-4	4ml
---------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Quantity Sufficient for 40 reactions of 0.1 ml each.

Technical Info / Product For the Manual please [click here](#).
Notes



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)