

Anti-Rabbit-AP-Polymer

For use in chromogenic in situ hybridization (CISH) procedures

The Anti-Rabbit-AP-Polymer (AB12) is intended to be used for detection steps in chromogenic in situ hybridization (CISH) applications on formalin-fixed, paraffin-embedded specimens. The Anti-Rabbit-AP-Polymer is intended to be used in combination with Digoxigenin-labeled ZytoFast CISH Probes and ZytoFast PLUS CISH Implementation Kits (T-1151-40 or T-1061-40).

Ordering Information

[Order Online »](#)

ZTV-AB-0012-4	4ml
---------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Species Reactivity Rabbit

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