

Anti-Mouse-HRP- Polymer

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

Polymerized HRP-conjugated secondary antibodies for the detection of mouse primary antibodies in chromogenic *in situ* hybridization (CISH). Component of *ZytoDot* 2C CISH Polymer Detection Kit (Prod. No. [ZTV-C-3028](#)).

Ordering Information

[Order Online »](#)

ZTV-AB-0002-4	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 Antibody to Horseradish peroxidase

Formulation Solution of salt (containing synthetic polypeptides) in water

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