

Nuclear Red Solution

For use in chromogenic in situ hybridization (CISH) procedures

The Nuclear Red Solution (CS3) is intended to be used for counterstaining steps in chromogenic in situ hybridization (CISH) applications on formalin-fixed, paraffin-embedded specimens. The Nuclear Red Solution is intended to be used in combination with Digoxigenin-labeled ZytoFast Probes and the ZytoFast PLUS CISH Implementation Kit AP-NBT/BCIP (T-1061-40).

Ordering Information

[Order Online »](#)

ZTV-CS-0003-20	20ml
----------------	------

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 0.5 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)