

ZytoDot® HRP-Green Solution Set

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

Ordering Information

[Order Online »](#)

| | |
|----------------|-----------|
| ZTV-C-3039-100 | 100 tests |
|----------------|-----------|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Contents HRP-Green Solution A, 0.6 ml
HRP-Green Solution B, 10 ml
HRP-Green Solution C, 0.6 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