

ZytoLight® FISH- Tissue Implementation Kit

For fluorescence *in situ* hybridization (FISH)
using any *ZytoLight* FISH probe

Ordering Information

[Order Online »](#)

ZTV-Z-2028-20	20 tests
---------------	----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Protect from light.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Contents Heat Pretreatment Solution Citric; Pepsin Solution; Wash Buffer SSC; 25x Wash Buffer A; DAPI/Antifade-Solution, good for 5 or 20 tests.

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