

25X Wash Buffer A

For use in fluorescence in situ hybridization
(FISH) procedures

Ordering Information

[Order Online »](#)

ZTV-WB-0002-50	50ml
----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Store, tightly closed, in the original vessel at temperatures between 2°C and 8°C
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance	liquid; colorless
Formulation	Solution of salt in water
Quantity	Sufficient for 17 staining jars of 70 ml each.
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