

# Cytology Pepsin Solution

For use in fluorescence in situ hybridization procedures on cytology specimens

## Ordering Information

[Order Online »](#)

ZTV-ES-0002-50	50ml
----------------	------

## Manuals, SDS & CofA

[View Online »](#)

FISH Accessories

## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Appearance</b>	liquid; colorless
<b>Formulation</b>	Solution of salt in water
<b>Quantity</b>	Sufficient for 40 reactions

**Technical Info / Product Notes** For the Manual please [click here](#).



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)