



# AMPIVIEW<sup>®</sup>

## Hybridization Buffer (1X)

### Optimized buffer to use with Enzo's AMPIVIEW<sup>®</sup> RNA probes

AMPIVIEW<sup>®</sup> Hybridization Buffer (1X) is a dilution or conditioning buffer used to dilute AMPIVIEW<sup>®</sup> RNA probes or precondition tissue samples prior to hybridization.

### Ordering Information

[Order Online »](#)

ENZ-ACC152-0010	10ml
-----------------	------

### Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Use/Stability** Stable for at least one year after receipt when stored as recommended. Following dilution in appropriate buffer system, product is stable for at least 2 days.

**Handling** Allow contents to warm up to room temperature prior to use.

**Short Term Storage** +4°C

**Long Term Storage** -20°C

**Shipping** Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

**Technical Info / Product Notes** Diluent buffer used for optimizing Enzo's AMPIVIEW<sup>®</sup> RNA Probes for ISH application.



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)