



# AMPIVIEW™ Hybridization Buffer (1X)

**REF** ENZ-ACC152-0010

10.0 mL AMPIVIEW™ Hybridization Buffer (1X)

## INTENDED USE:

For Research Use Only. Not for use in diagnostic procedures.

## SUMMARY AND EXPLANATION

AMPIVIEW™ Hybridization Buffer (1X) is a ready-to-use formamide-based buffer optimized for use with AMPIVIEW™ RNA probes, accessory and detection reagents to achieve appropriate *in situ* hybridization (ISH) staining. During ISH process, AMPIVIEW Hybridization Buffer (1X) is used to dilute probes and create appropriate hybridization conditions on the slide.

## KNOWN APPLICATION

*in situ* Hybridization (ISH) on formalin-fixed paraffin-embedded (FFPE) tissue specimens or cells.

## PRODUCTS SUPPLIED

AMPIVIEW™ Hybridization Buffer (1X) (ENZ-ACC152)

## STORAGE AND SHELF-LIFE

Upon receipt, store probes solution and hybridization buffer at -20°C. This product is stable under these conditions up to the expiration date indicated in the vial label.

## PERFORMANCE CONSIDERATIONS

1. Do not use reagents past their expiration date.
2. Do not allow the slides to dry completely during the hybridization and detection procedures, or erroneous results may occur. Avoid drying by ensuring that the entire specimen is covered with sufficient amounts of buffers and reagents as recommended in the procedures. While incubating, the slides may be covered with a cover slip to help prevent drying.
3. Cross-contamination of samples could cause false results. Use care preparing slides for more than one specimen.
4. Allow all components to reach room temperature (20-30°C) before beginning the test procedure.
5. Incubation times and temperatures other than those specified may give erroneous results.
6. Improper specimen preparation may cause false results.

## LIMITATIONS

This procedure is for research use only. It is not intended for diagnostic or therapeutic use.

## PRECAUTIONS

1. Refer to reagent Safety Data Sheet (SDS) for precautions.
2. Specimens, before and after fixation, and all materials exposed to them should be handled and disposed of with proper precautions.
3. Never pipette reagents by mouth and avoid contact with skin and mucous membranes with reagents and specimens. If reagents and/or specimens come into contact with sensitive areas, rinse thoroughly with water and follow your institution's safety protocols.
4. Avoid RNases, which are commonly found on skin. Wear gloves at all times, and work in a clean environment.

## TECHNICAL NOTES

1. AMPIVIEW™ Hybridization buffer is a ready-to-use dilution or conditioning buffer.
2. AMPIVIEW™ RNA probes are diluted in AMPIVIEW™ Hybridization buffer.
3. AMPIVIEW™ RNA probes have been optimized with DIGX® anti-digoxigenin (mouse or rabbit) and POLYVIEW® PLUS Detection Systems in combination with HIGHDEF® chromogens. Use of other linkers and polymer-based reagents will require prior optimization.
4. Autostainers such as LEICA BOND, Roche DISCOVERY or other automated systems can be used, but will require prior optimization. AMPIVIEW™ RNA probes might need to be diluted with the hybridization buffer for automated systems.
5. Changes in the amount of probe or temperature incubation times from what is recommended without prior testing, may lead to inconsistent results.

For Research Use Only, Not for Use in Diagnostics Procedures.

All product names, logos, brands and trademarks are registered and legally protected by their respective owners.

PM-PL0099 Rev101923

### GLOBAL HEADQUARTERS

Enzo Life Sciences, Inc.  
81 Executive Blvd, Ste 3  
Farmingdale, NY 11735 USA  
T 1-800-942-0430  
F 1-610-941-9252  
E info-usa@enzolifesciences.com  
www.enzolifesciences.com

### EUROPE

Enzo Life Sciences (ELS) AG  
Industriestrasse 17, Postfach  
CH-4415 Lausen, Switzerland  
T +41 61 926 89 89  
F +41 61 926 89 79  
E info-ch@enzolifesciences.com  
www.enzolifesciences.com