

# AMPIVIEW<sup>®</sup> Wash Buffers Kit

Optimized wash buffer 1 and wash buffer 2 to use with Enzo's AMPIVIEW<sup>®</sup> RNA probes

AMPIVIEW<sup>®</sup> Wash Buffers Kit contains AMPIVIEW<sup>®</sup> Wash Buffer 1 and AMPIVIEW<sup>®</sup> Wash Buffer 2, which are used for post-hybridization washes prior to detection. AMPIVIEW<sup>®</sup> Wash Buffer 1 is used to remove all non-bound AMPIVIEW<sup>®</sup> RNA Probes and AMPIVIEW<sup>®</sup> Wash Buffer 2 is used to remove Wash Buffer 1 before the detection step.

## Ordering Information

[Order Online »](#)

ENZ-ACC161-0040	40ml
-----------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	Stable for at least one year after receipt when stored as recommended.
Handling	Allow contents to warm up to 37°C prior to use.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Contents	1 x 20 mL AMPIVIEW <sup>®</sup> Wash Buffer 1 1 x 20 mL AMPIVIEW <sup>®</sup> Wash Buffer 2
----------	--

Technical Info / Product Notes	Wash buffers to remove non-bound probes to prevent background or non-specific detection.
--------------------------------	--



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)