

AMPIVIEW® Wash Buffers Kit

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

AMPIVIEW® Wash Buffers Kit contains AMPIVIEW® Wash Buffer 1 and AMPIVIEW® Wash Buffer 2, which are used for post-hybridization washes prior to detection. AMPIVIEW® Wash Buffer 1 is used to remove all non-bound AMPIVIEW® RNA Probes and AMPIVIEW® 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® Wash Buffer 1
1 x 20 mL AMPIVIEW® Wash Buffer 2

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

Last modified: October 17, 2024



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