

AMPIVIEW® Wash Buffer (for Ambient Use)

Ready-to-use post-hybridization wash buffer for room-temperature ISH workflows; A non-formamide solution for manual and automated systems that would require ambient workflow

AMPIVIEW® Wash Buffer (for Ambient Use) is a ready-to-use post-hybridization wash solution for AMPIVIEW® RNA in situ hybridization (ISH) workflows performed at room temperature, complementing the heated wash steps used with the standard AMPIVIEW® Wash Buffer Kit (ENZ-ACC161).

Ordering Information

[Order Online »](#)

ENZ-ACC188-0250	250ml
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

- Ready-to-use, no preparation required
- Formulated for ambient-temperature (room-temperature) wash steps
- Compatible with AMPIVIEW® RNA probe ISH workflows

Handling & Storage

Use/Stability	Stable under recommended storage conditions up to the expiration date indicated on the vial label.
Handling	Allow to reach room temperature before use.
Short Term Storage	Room Temperature
Long Term Storage	2–8°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Contents	AMPIVIEW® Wash Buffer (for Ambient Use) (ENZ-ACC188)
Technical Info / Product Notes	Used during post-hybridization wash steps in the AMPIVIEW® ISH protocol, as an ambient-temperature alternative to the heated (45°C) wash buffer steps. Note: no product insert/IFU was available for this SKU at time of import; content drafted from AMPIVIEW® product-line conventions and should be verified against the official insert before publishing.



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)