

AMPIVIEW®

Hybridization Buffer (Fast Hyb Solution)

Ready-to-use hybridization buffer for AMPIVIEW® RNA probes; Fast hybridization solution

AMPIVIEW® Hybridization Buffer (Fast Hyb Solution) is a ready-to-use buffer optimized for use with AMPIVIEW® probes, accessory, and detection reagents to achieve appropriate in situ hybridization (ISH) staining. During the ISH process, this buffer is used to dilute probes and create Optimal hybridization conditions .

Supplied as 1 x 100 mL.

Ordering Information

[Order Online »](#)

ENZ-ACC187-0100	100ml
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

- Ready-to-use dilution/conditioning buffer — no preparation required
- Formamide-based formula optimized for AMPIVIEW® RNA probes
- Compatible with manual and automated ISH staining platforms (prior optimization required for autostainers)

Handling & Storage

Use/Stability	Stable under recommended storage conditions up to the expiration date indicated on the vial label.
Handling	Do not allow slides to dry during hybridization/detection. Allow components to reach room temperature (20–30°C) before use.
Short Term Storage	2–8°C
Long Term Storage	2–8°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Contents	1 x 100 mL AMPIVIEW® Hybridization Buffer (1X, Fast Hyb Solution)
Technical Info / Product Notes	AMPIVIEW® RNA probes are diluted in AMPIVIEW® Hybridization Buffer — a fast hybridization solution optimized to promote molecular crowding, reduce background, and achieve highly specific probe binding



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)