

AMPIVIEW® Hybridization Buffer (1X)

Optimized buffer to use with Enzo's
AMPIVIEW® RNA probes

AMPIVIEW® Hybridization Buffer (1X) is a dilution or conditioning buffer used to dilute AMPIVIEW® RNA probes or precondition tissue samples prior to hybridization.

Ordering Information

[Order Online »](#)

ENZ-ACC152-0010	10ml
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability Stable for at least one year after receipt when stored as recommended. Following dilution in appropriate buffer system, product is stable for at least 2 days.

Handling Allow contents to warm up to room temperature prior to use.

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Technical Info / Product Notes Diluent buffer used for optimizing Enzo's AMPIVIEW[®] RNA Probes for ISH application.

Last modified: October 17, 2024



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)