

In situ hybridization buffer (1.25X concentrate)

In Situ Hybridization Buffer provides a high quality hybridization medium for *in situ* hybridization analyses using either DNA or oligonucleotide probes labeled with biotin-, digoxigenin- or fluorescein-modified nucleotides.

Citations: 14

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-33808	10ml
-----------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability Stable for up to 1 year when stored refrigerated in a well sealed bottle.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Formulation Liquid. Sodium phosphate, NaCl and EDTA (SSPE) buffered formamide, containing hybridization enhancers.

Quantity Provides a sufficient amount of hybridization buffer concentrate to prepare 12.5ml of ready-to-use probe reagent solutions.

Last modified: May 29, 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)