

HighFidelity LNA PCR Labeling Kit

HighFidelity LNA PCR Labeling Kit is designed to produce randomly Locked nucleic acid (LNA)-modified DNA probes by PCR. Such probes are ideally suited for in situ hybridization and Northern Blot experiments. The ribose ring of locked nucleic acids is "locked" in the ideal conformation for Watson-Crick binding. LNA-modified DNA probes therefore possess an unprecedented thermal stability upon hybridization resulting in an increased sensitivity. LNA-UTP, LNA-CTP, LNA-ATP and LNA-GTP are efficiently incorporated into DNA as substitute for their natural counterpart (dTTP, dCTP, dATP or dGTP, respectively) using an optimized reaction buffer and a High Fidelity Polymerase. 50 % LNA-NTP substitution typically results in an optimal balance between reaction and labeling efficiency. Individual optimization of LNA-NTP/dNTP ratio however, can easily be achieved with the single nucleotide format. The kit contains sufficient reagents for 10 labeling reactions each LNA-NTP (20 µl each (50 % LNA-NTP substitution))."

Ordering Information

[Order Online »](#)

JBS-APP-101-LNA	40Reactions
-----------------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes Preparation of LNA-labeled DNA probes by PCR

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



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