

AMPINEXT™ NGS

Barcode (Index) Set-12

Designed to construct multiplex DNA/RNA libraries used for next-generation sequencing with the Illumina platform

The AMPINEXT™ NGS Barcode (Index) Set-12 is designed to construct multiplex DNA/RNA libraries used for next-generation sequencing with the Illumina platform including GAllx, HiSeq, MiSeq, and NextSeq. This product includes 12 DNA barcodes, each of which contain an index and flow cell binding sequence and can be attached to the sample insert during library preparation. Multiplexing several samples in a single flow cell can be achieved by pooling with the AMPINEXT™ NGS Barcodes. Pooling barcoded samples into a single flow cell significantly reduces hands-on time and provides robust data quality in NGS.

Manuals, SDS & CofA

[View Online »](#)

- Compatible with various library preparation workflows.
- High throughput and cost reduction: Up to 144 multiplexed samples can be processed.
- The indices are designed to be error resistant and allow for proper differentiation between samples.
- Minimizes PCR amplification-based bias and prevents poor reads from single base errors introduced during PCR.

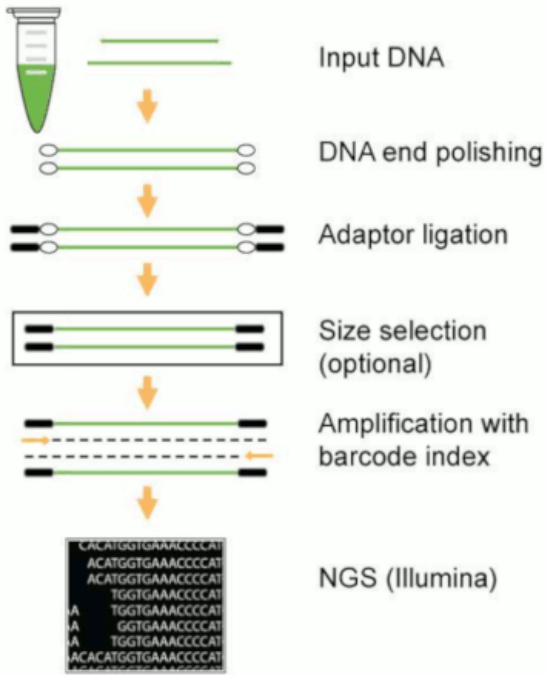


Figure 1. Schematic Procedure for using the AMPINEXT NGS Barcode (Index) Set-12

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes

Designed to construct multiplex DNA/RNA libraries used for next-generation sequencing with the Illumina platform

Contents

AMPINEXT™ Barcodes 1-12 (10 µM)
AMPINEXT™ Universal Primer (10 µM)

Last modified: May 29, 2024



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