

## Tools to Streamline Your NGS Workflow

Next generation sequencing (NGS) is a method of high-throughput sequencing in which millions of DNA strands can be sequenced in parallel. Many researchers use NGS as an important tool to gain insight into the detailed and comprehensive genetic and epigenetic modifications of genomes from many different cell types. Choose from our comprehensive mix of kits to extract and prepare your library sample for NGS analysis. Whether you are preparing samples for ChIP-seq, whole genome sequencing, or Bisulfite-sequencing our kits are easy, efficient, and convenient.

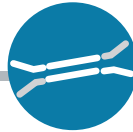
### Reliable and Streamlined NGS Workflow

#### Nucleic Acid Extraction



AMPIXTRACT® Solutions

#### Library Preparation



AMPINEXT™ Products

#### Sequencing & Analysis



Compatible with Illumina® Platform

#### Nucleic Acid Extraction



- **Fast** – Entire procedure can be finished in as little as 15 minutes with consistent isolation conditions
- **Effective** – High efficiency of nucleic acid isolation from various sample types with low amounts of start material (as low as 1 ng)
- **Reliable** – Our specially designed columns allow for nucleic acids to be conveniently recovered
- **Safe** – Use of non-toxic reagents and no phenol chloroform

#### Library Preparation

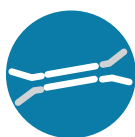


- **Fast and Streamlined Procedure** – Compact protocols with simple clean-up steps to save time and prevent handling errors, as well as loss of valuable samples
- **Highly Convenient** – The kit contains all required components for each step of the library preparation for the most reliable and consistent results
- **Minimize bias** – High quality enzymes achieve reproducibly high yields of nucleic acid library with minimal sequence bias and low error rates
- **Flexibility** – Can be used for both non-barcoded (singleplexed) and barcoded (multiplexed) library preparation with broad range of nucleic acid input



## NUCLEIC ACID EXTRACTION

Product Name	Product #	Sample Input	Elution Volume	Protocol Time
AMPIXTRACT® General Tissue Section DNA Isolation Kit	ENZ-GEN500	1 ng (optimal range 10 ng – 1 µg)	8 - 20 µL	2 hours
AMPIXTRACT® Paraffin Tissue Section DNA Isolation Kit	ENZ-GEN501	1 ng (optimal range 10 ng – 1 µg)	8 - 20 µL	2 hours
AMPIXTRACT® Urine DNA Isolation Kit	ENZ-GEN502	1 ng – 2 µg (optimal range 10 ng – 1 µg)	8 - 20 µL	20 minutes
AMPIXTRACT® Blood and Cultured Cell DNA Extraction Kit	ENZ-GEN503	1 ng – 4 µg (optimal range 10 ng – 1 µg)	10 - 20 µL	20 minutes
AMPIXTRACT® DNA Isolation Kit for Plasma/Serum	ENZ-45018	0.1 ng (optimal range 1 ng – 1 µg)	8 - 20 µL	15 minutes



## LIBRARY PREPARATION

Product Name	DNA		Bisulfite-Seq			Post Bisulfite-Seq	ChIP-Seq	
		AMPINEXT™ DNA Library Preparation Kit	AMPINEXT™ High-Sensitivity DNA Library Preparation Kit	AMPINEXT™ High-Sensitivity Bisulfite-Seq Kit	AMPINEXT™ RNA Bisulfite-Seq Kit	AMPINEXT™ 5-mC RNA Bisulfite-Seq Easy Kit	AMPINEXT™ Post-Bisulfite DNA Library Preparation Kit	AMPINEXT™ ChIP-Seq High-Sensitivity Kit
Product #	ENZ-GEN504	ENZ-GEN505	ENZ-GEN508	ENZ-GEN510	ENZ-GEN511	ENZ-GEN512	ENZ-GEN509	
Input Type	DNA	DNA	DNA	RNA	RNA	DNA	Cells	Tissues
Input Range (Optimal Range)	5 ng – 1 µg (100 ng – 200 ng)	0.2 ng – 100 ng (10 ng – 50 ng)	0.2 ng – 200 ng (10 ng – 50 ng)	5 ng – 1 µg (200 ng – 500 ng)	5 ng – 500 ng (100 ng – 200 ng)	0.5 ng – 1 µg (100 ng – 200 ng)	10 <sup>5</sup> – 10 <sup>6</sup> Cells  (4×10 <sup>5</sup> – 5×10 <sup>5</sup> ) Cells	5 mg – 50 mg  (20 mg – 30 mg)
Protocol Time	< 2.5 hours	< 1.5 hours	< 6.5 hours	6 hours		5 hours	< 7 hours	

## RELATED PRODUCTS

Product Name	Product Number	Contents
AMPINEXT™ DNA Size Selection Kit	ENZ-GEN506	MQ Binding Beads, Elution Solution, 96-well plate
AMPINEXT™ NGS Barcode (Index) Set-12	ENZ-GEN507	AMPINEXT™ Barcodes 1-12 (10 µM) AMPINEXT™ Universal Primer (10 µM)

All product names, logos, and brands are the property of their respective owners.

**Global Headquarters**  
**ENZO LIFE SCIENCES, INC.**  
 Ph: 800.942.0430  
 info-usa@enzolifesciences.com

**European Sales Office**  
**ENZO LIFE SCIENCES (ELS) AG**  
 Ph: +41 61 926 8989  
 info-eu@enzolifesciences.com

**Belgium, The Netherlands & Luxembourg**  
**& Luxembourg**  
 Ph: +32 3 466 0420  
 info-be@enzolifesciences.com

**France**  
 Ph: +33 472 440 655  
 info-fr@enzolifesciences.com

**Germany**  
 Ph: +49 7621 5500 526  
 info-de@enzolifesciences.com

**UK & Ireland**  
 Ph: 0845 601 1488 (UK customers)  
 Ph: +44 1392 825900  
 info-uk@enzolifesciences.com