

GenNature Adapter Set Quick Manual

(Cat. No.: NGS01204)

For Illumina Sequencing Platform

Adapter information

GN701–GN707 Adapters

CAAGCAGAAGACGGCATACGAGAT[i7]GTCTCGTGGGCTCGG

GS505–GS506 Adapters

AATGATACGGCGACCACCGAGATCTACAC[i5]TCGTCGGCAGCGTC

GN	Bases in GN	i7 Bases for Sample Sheet	GS	Bases in GS	i5 Bases for Sample Sheet *	i5 Bases for Sample Sheet **
GN701	TCGCCTTA	TAAGGCGA	GS505	GTAAGGAG	GTAAGGAG	CTCCTTAC
GN702	CTAGTACG	CGTACTAG	GS506	ACTGCATA	ACTGCATA	TATGCAGT
GN704	GCTCAGGA	TCCTGAGC				
GN705	AGGAGTCC	GGACTCCT				
GN706	CATGCCTA	TAGGCATG				
GN707	GTAGAGAG	CTCTCTAC				

* Compatible with Illumina sequencer, NovaSeq, MiSeq, and HiSeq 2000/2500

** Compatible with Illumina sequencer, iSeq, MiniSeq, NextSeq, and HiSeq 3000/4000

Remain stable for 12 months. Avoid repeated freeze-thaw cycles.

Pooling Guidelines

Single Indexing

The following table shows strategies for creating 2~12-plex pools of dual-indexed libraries for single-indexed sequencing, which sequences Index 1 only.

Plexity	GN	GS
2	• GN701 and GN702	Any GS
3	• GN701, GN702, and GN704	Any GS
4 or 5	• GN701, GN702, GN704 and any other GN	Any GS
6-12	• GN701, GN702, GN704, GN705, GN706 and GN707	Any GS

Pooling Guidelines

Dual Indexing

The following table shows strategies for pooling dual-indexed libraries for dual-indexed sequencing, which sequences Index 1 and Index 2.

Plexity	GN	GS
2–6	At least two unique GN	Two unique GS
7–12	One of the following combinations: <ul style="list-style-type: none">• GN701, GN702, GN704 and any other GN• GN705, GN706, GN707 and any other GN	GS505 and GS506

Welcome to download detailed user guide from [GenScript official website](#).

Version 2019-08-1