

Catalog No.: 53155

**Description:** Kit contains 100 µL of each primer including 8 i7 Indexing Primers and 12 i5 Indexing Primers for multiplexing up to 96 samples for sequencing on Illumina Next-Generation Sequencing Systems. Oligonucleotide sequences ©2019 Illumina Inc. All rights reserved.

**Store at:** -20°C

Primers are supplied as 25 µM in 0.1 mM low EDTA TE Buffer for use in multiplexing samples generated in kits such as the ATAC-Seq Kit (Cat. No. 53150) and CUT&Tag-IT™ Assay Kits (Cat. Nos. 53160, 53165)

## Index Primers and Sample Sheet Information

Index 1 (i7) Primers

CAAGCAGAAGACGGCATAACGAGAT[(i7)GTCTCGTGGGCTCGG

Index 2 (i5) Primers

AATGATACGGCGACCACCGAGATCTACAC[i5]TCGTCCGGCAGCGTC

i7 Index Primer Name	i7 Index Name	i7 Sequence	Sample sheet
i7 Indexed Primer 1	N701	TCGCCTTA	TAAGGCCGA
i7 Indexed Primer 2	N702	CTAGTACG	CGTACTAG
i7 Indexed Primer 3	N703	TTCTGCCT	AGGCAGAA
i7 Indexed Primer 4	N704	GCTCAGGA	TCCTGAGC
i7 Indexed Primer 5	N705	AGGAGTCC	GGACTCCT
i7 Indexed Primer 6	N706	CATGCCTA	TAGGCATG
i7 Indexed Primer 7	N707	GTAGAGAG	CTCTCTAC
i7 Indexed Primer 8	N708	CCTCTCTG	CAGAGAGG

**Sample Sheet Information for NovaSeq v 1.0 Reagent Kits, MiSeq, and HiSeq 2000/2500)**

i5 Index Primer Name	i5 Index Name	i5 Sequence	Sample sheet
i5 Indexed Primer 1	N501	TAGATCGC	TAGATCGC
i5 Indexed Primer 2	N502	CTCTCTAT	CTCTCTAT
i5 Indexed Primer 3	N503	TATCCTCT	TATCCTCT
i5 Indexed Primer 4	N504	AGAGTAGA	AGAGTAGA
i5 Indexed Primer 5	N505	GTAAGGAG	GTAAGGAG
i5 Indexed Primer 6	N506	ACTGCATA	ACTGCATA
i5 Indexed Primer 7	N507	AAGGAGTA	AAGGAGTA
i5 Indexed Primer 8	N508	CTAAGCCT	CTAAGCCT
i5 Indexed Primer 10	N510	CGTCTAAT	CGTCTAAT
i5 Indexed Primer 11	N511	TCTCTCCG	TCTCTCCG
i5 Indexed Primer 13	N513	TCGACTAG	TCGACTAG
i5 Indexed Primer 15	N515	TTCTAGCT	TTCTAGCT

**Sample Sheet Information for NovaSeq v 1.5 Reagent Kits, iSeq, MiniSeq, NextSeq, and HiSeq 3000/4000)**

i5 Index Primer Name	i5 Index Name	i5 Sequence	Sample sheet
i5 Indexed Primer 1	N501	TAGATCGC	GCGATCTA
i5 Indexed Primer 2	N502	CTCTCTAT	ATAGAGAG
i5 Indexed Primer 3	N503	TATCCTCT	AGAGGATA
i5 Indexed Primer 4	N504	AGAGTAGA	TCTACTCT
i5 Indexed Primer 5	N505	GTAAGGAG	CTCCTTAC
i5 Indexed Primer 6	N506	ACTGCATA	TATGCAGT
i5 Indexed Primer 7	N507	AAGGAGTA	TACTCCTT
i5 Indexed Primer 8	N508	CTAAGCCT	AGGCTTAG
i5 Indexed Primer 10	N510	CGTCTAAT	ATTAGACG
i5 Indexed Primer 11	N511	TCTCTCCG	CGGAGAGA
i5 Indexed Primer 13	N513	TCGACTAG	CTAGTCGA
i5 Indexed Primer 15	N515	TTCTAGCT	AGCTAGAA