

## Dual Index Primers Set 1 for Illumina®

---

**Catalog No.:** 53221

**Description:** Kit contains 10  $\mu$ M stocks of each primer including 20  $\mu$ L each of 6 i5 Indexing Primers and 15  $\mu$ L each of 8 i7 Indexing Primers for multiplexing up to 48 samples generated with DNA Library Prep Kit for Illumina® (Catalog No. 53220) for sequencing on Illumina Next-Generation Sequencing Systems. Can be combined with Dual Index Primers Set 2 for Illumina® (Catalog No. 53222) to multiplex up to 96 samples. Full primer sequences can be downloaded from the product page at [activemotif.com](http://activemotif.com). Oligonucleotide sequences ©2019 Illumina Inc. All rights reserved.

**Store at:** -20°C

### Kit Contents:

Index Primer Name	Amount
AM i5-001 Primer	20 $\mu$ L
AM i5-002 Primer	20 $\mu$ L
AM i5-003 Primer	20 $\mu$ L
AM i5-004 Primer	20 $\mu$ L
AM i5-005 Primer	20 $\mu$ L
AM i5-006 Primer	20 $\mu$ L
AM i7-001 Primer	15 $\mu$ L
AM i7-002 Primer	15 $\mu$ L
AM i7-003 Primer	15 $\mu$ L
AM i7-004 Primer	15 $\mu$ L
AM i7-005 Primer	15 $\mu$ L
AM i7-006 Primer	15 $\mu$ L
AM i7-007 Primer	15 $\mu$ L
AM i7-008 Primer	15 $\mu$ L

## Index Primers and Sample Sheet Information

### i5 Primer

5'-AATGATACGGCGACCACCGAGATCTACAC[i5]ACACTCTTCCCTACACGACGCTCTTCCGATC\*T-3'

### i7 Primer

5'-CAAGCAGAAGACGGCATAACGAGAT[i7]GTGACTGGAGTTCAGACGTGTGCTCTTCCGA\*T-3'

### Sample Sheet Information

Index Primer Name	Sample sheet for MiSeq, HiSeq 200/2500, Nova-Seq v1.0 Reagent Kits	Sample sheet for iSeq, MiniSeq, NextSeq, HiSeq 3000/4000, Nova-Seq v1.5 Reagent Kits
<b>i5 Primers</b>		
AM i5-001 Primer	AGTTGAAT	ATTCAACT
AM i5-002 Primer	ACCGGCCA	TGGCCGGT
AM i5-003 Primer	CTGAACCG	TATCCTCT
AM i5-004 Primer	GCGTGCTC	GAGCACGC
AM i5-005 Primer	ATACCGTT	AACGGTAT
AM i5-006 Primer	CGGTCCTA	TAGGACCG
<b>i7 Primers</b>		
AM i7-001 Primer	GCAGTCTT	GCAGTCTT
AM i7-002 Primer	TGATCAGT	TGATCAGT
AM i7-003 Primer	GTCGGCAC	GTCGGCAC
AM i7-004 Primer	TGGAAGAG	TGGAAGAG
AM i7-005 Primer	CGGAAGGT	CGGAAGGT
AM i7-006 Primer	AATCTGGT	AATCTGGT
AM i7-007 Primer	ATGTGCCT	ATGTGCCT
AM i7-008 Primer	GTGTCCTG	GTGTCCTG