

Nextera™-Compatible UDI Primers (1 - 96)

Catalog No: 53178

Format: 1 kit

Background: Kit contains 10 µL of pre-mixed unique dual indexed primers for multiplexing up to 96 samples for sequencing on Illumina Next-Generation Sequencing Systems. Oligonucleotide sequences ©2019 Illumina Inc. All rights reserved.

Primers are supplied as 12.5 µM each (25 µM combined) 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)TCGTCGGCAGCGTC

For the full plate map with index sequences, please visit the product web page at www.activemotif.com.

Quality Control: Unique Dual Index (UDI) primers are subject to rigorous quality control testing, including sequence verification, purity assessment, concentration confirmation, and functional performance validation, to ensure accurate sample indexing and consistent sequencing results.

Storage: Store at -20°C.

Guarantee: This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.