

CUTANA™ Nextera-Compatible CDI Primer Set 1

Catalog No	14-1190-24rxn	Pack Size	24 Reactions
Lot No	26174001-52	Kit Version	v1

DESCRIPTION

CUTANA™ Nextera-Compatible CDI Primer Set 1 provides purpose-built, combinatorial dual-index (CDI) primers for generating indexed, sequencing-ready libraries from ATAC-seq and CUT&Tag workflows. The primers are supplied ready to drop into CUTANA™ ATAC-seq and CUTANA™ CUT&Tag assays with no protocol changes. The set contains 4x i5 and 6x i7 indices, yielding 24 combinatorial dual-index libraries per set.

Key features

- Purpose-built dual-index primers for ATAC-seq and CUT&Tag
- Ready to use out of the box: no index design, ordering, or titration required
- Drops into CUTANA™ ATAC-seq and CUT&Tag with no protocol changes
- Fully compatible with Tn5-generated libraries and standard Illumina sequencing

For the complete ATAC-seq workflow, pair with the CUTANA™ ATAC-seq Kit (14-1121) or the CUTANA™ Tn5 Transposome (15-1036).

SET CONTENTS

<u>Item</u>	<u>Cat No</u>	<u>Item</u>	<u>Cat No</u>
i701 Primer	18-1301-13	i501 Primer	18-1201-13
i702 Primer	18-1302-13	i502 Primer	18-1202-13
i703 Primer	18-1303-13	i503 Primer	18-1203-13
i704 Primer	18-1304-13	i504 Primer	18-1204-13
i705 Primer	18-1305-13		
i706 Primer	18-1306-13		

RECOMMENDED ACCESSORY PRODUCTS

ATAC-seq | Kit-based workflow

<u>Item</u>	<u>Cat No</u>
CUTANA™ ATAC-seq Kit	14-1121

ATAC-seq | User-defined Tn5-based workflow

<u>Item</u>	<u>Cat No</u>
CUTANA™ Tn5 Transposome	15-1036
Non-Hot Start 2X PCR Master Mix	15-1018

CUT&Tag | User-defined pAG-Tn5 workflow

<u>Item</u>	<u>Cat No</u>
CUTANA™ pAG Tn5 for CUT&Tag	15-1017
Non-Hot Start 2X PCR Master Mix	15-1018

TECHNICAL INFORMATION

Storage	Store at -20°C. Refer to product label for expiry date.
Instructions for Use	Each CUTANA™ Nextera-Compatible CDI Primer Set 1 produces 24 combinatorial dual-index libraries, with enough primer volume for 1x 24 ATAC-seq reactions or 3x 24 CUT&Tag reactions.

CERTIFICATE OF ANALYSIS

Quality Control Metrics

Result

Concentration and Purity (NanoDrop)

Pass

Each primer in the set is individually quantified by UV absorbance (NanoDrop) to confirm concentration, with purity assessed by 260/280 and 260/230 absorbance ratios.

Sequence Identity (Mass Spectrometry)

Pass

Each primer in the set is individually analyzed by mass spectrometry to confirm sequence identity (observed mass matches expected mass).

Barcode Identity and Cross-Contamination (Cloning and Sequencing)

Pass

Each lot is validated for barcode identity and absence of cross-contamination through amplification of a defined template, cloning, and sequencing to confirm the expected barcode identity of each primer and the absence of cross-contamination between primers.

Functional Testing (ATAC-seq Library Construction)

Pass

Each lot is functionally validated by amplification of ATAC-seq libraries generated from 50,000 K562 nuclei alongside established EpiCypher references. Libraries are sequenced on an Illumina NextSeq2000 with paired-end sequencing (2x50 bp). Assessment is based on metrics including library yields and fragment size distribution, percent unique reads, percent mitochondrial reads, percent *E. coli* reads, fraction of reads in peaks (FRiP), insert size distribution, and TSS enrichment score.

PUBLISH DATE

9/15/2026

APPROVED BY



ANDREA JOHNSTONE, PHD

SENIOR DIRECTOR OF PRODUCT MANAGEMENT

This product ("Product") is warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications, and/or accompanying package inserts ("Documentation") and to be free from defects in material and workmanship. Unless otherwise expressly authorized in writing, Products are supplied for research use only. No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper, and intended usage. This warranty does not extend to anyone other than the original purchase of the Product ("Buyer"). No other warranties, express or implied, are granted, including without limitation, implied warranties of merchantability, fitness for any particular purpose, or non infringement. Buyer's exclusive remedy for non-conforming Products during the warranty period is limited to replacement of or refund for the non-conforming Product(s). There is no obligation to replace Products as the result of (i) accident, disaster or event of force majeure, (ii) misuse, fault or negligence of or by Buyer, (iii) use of the Products in a manner for which they were not designed, or (iv) improper storage and handling of the Products.