

CUTANA™ CUT&RUN Library Prep Kit with Primer Set 2

Catalog No	14-1001-48s2	Pack Size	48 Reactions
Lot No	26047007-51	Kit Version	v1

DESCRIPTION

The CUTANA™ CUT&RUN Library Prep Kit offers high fidelity library generation for Illumina® sequencing by harnessing the power of New England Biolabs® best-in-class NEBNext® reagents. The kit offers a streamlined protocol specifically optimized for high sensitivity CUT&RUN applications, including those with low cell inputs. Included are all necessary reagents to perform end repair, adaptor ligation, combinatorial dual indexing for multiplexing up to 48 samples, and DNA cleanup with CUTANA™ DNA Purification Beads (EpiCypher 21-1407). If additional multiplexing is desired, this kit can be used in tandem with Primer Set 1 (EpiCypher 14-1001-48s1) for up to 96 samples. Pairing this kit with the CUTANA™ ChIC/CUT&RUN Kit (EpiCypher 14-1048) affords users a cells-to-sequencing solution for chromatin mapping experiments with all the necessary controls and validated reagents to ensure confidence in obtaining high quality data.

KIT CONTENTS

<u>Item</u>	<u>Cat No</u>	<u>Item</u>	<u>Cat No</u>
8-strip Tubes	10-0009p	Adapter for Illumina®	18-1000p
CUTANA™ DNA Purification Beads	21-1407p	Ligation Mix	15-1020p
0.1X TE Buffer	21-1025p	Ligation Enhancer	15-1021p
End Prep Enzyme	15-1019p	U-Excision Enzyme	15-1023p
End Prep Buffer	21-1012p	Hot-Start 2X PCR Master Mix	15-1022p
Multiplexing Primers	Primer Set 2 includes i5 primers 1-8 and i7 primers 7-12		

TECHNICAL INFORMATION

Storage	OPEN KIT IMMEDIATELY and store components at room temperature and -20°C as indicated (see Kit Manual for full instructions). Stable for 6 months upon date of receipt.
Instructions for Use	See the included Quick-Start Card. Find the User Manual and Quick-Start Card online in the “Resources” section at epicypher.com/14-1001 .

CERTIFICATE OF ANALYSIS

Quality Control Metrics

Functional Testing (CUT&RUN)

Each lot of CUTANA™ CUT&RUN Library Prep Kit is functionally validated by comparison to a previous lot and to established EpiCypher internal controls. CUT&RUN is performed using the CUTANA™ ChIC/CUT&RUN Kit (EpiCypher 14-1048) on 500k native K562 cells with 0.5 µg of IgG (EpiCypher 13-0042), H3K4me3 (EpiCypher 13-0060), H3K27me3 (EpiCypher 13-0055), or 0.125 µg CTCF (EpiCypher 13-2014) antibodies. Library preparation using the CUTANA™ CUT&RUN Library Prep Kit is performed with 0.5 - 5 ng inputs of CUT&RUN prepared DNA to assess performance across a range of input amounts. Libraries are sequenced on an Illumina NextSeq2000 with paired-end sequencing (2x50 bp). Assessment is based on metrics including DNA yields, library fragment size distribution, sequencing statistics, and genome-wide enrichment patterns.

Pass

PUBLISH DATE

6/9/2026

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.