

CUTANA™ ATAC-seq Kit

Catalog No	14-1121-24rxn	Pack Size	24 Reactions
Lot No	26107001-51	Kit Version	v1

DESCRIPTION

The **CUTANA™ ATAC-seq Kit** provides validated reagents and a streamlined workflow for generating sequencing-ready ATAC-seq libraries from cells or isolated nuclei. ATAC-seq was first described by Buenrostro et al. as a rapid method for mapping accessible chromatin by direct *in vitro* transposition of sequencing adapters into native chromatin [1]. The assay enables genome-wide analysis of open chromatin and regulatory element activity from low nuclei inputs.

The CUTANA™ ATAC-seq Kit includes pre-charged CUTANA™ Tn5 Transposome for ATAC-seq and optimized buffers for robust, reproducible chromatin accessibility profiling. Built on the foundation of the Omni-ATAC protocol described by Corces et al. and detailed by Grandi et al. [2,3], which established the performance standards researchers have come to expect from ATAC-seq, the CUTANA™ ATAC-seq Kit reimagines the workflow with optimized reagents and streamlined assay conditions designed to deliver a fully validated, user-friendly experience. These improvements support higher library yields, improved signal-to-noise, and consistent generation of high-quality sequencing-ready libraries from cells or isolated nuclei.

Key features

- Streamlined workflow from cells to sequencing-ready ATAC-seq libraries
- Includes pre-charged CUTANA™ Tn5 Transposome for ATAC-seq
- Built on proven Omni-ATAC workflow principles
- Optimized buffers and workflow conditions for robust, reproducible performance

For customers who prefer an end-to-end experience, see CUTANA™ ATAC-seq Services.

KIT CONTENTS

<u>Item</u>	<u>Cat No</u>	<u>Item</u>	<u>Cat No</u>
8-strip tubes	10-0009-12	Lysis Buffer	21-1039-12
10X PBS	21-1035-12	2X Tagmentation Buffer	21-1040-12
Stop Solution	21-1036-12	Enzyme Diluent	21-1041-12
DNA Purification Beads	21-1407-12	Non-Hot Start 2X PCR Master Mix	15-1018-12
0.1X TE Buffer	21-1025-12	5% Digitonin	21-1004-12
10X Wash Buffer	21-1037-12	Tn5 Transposome	15-1036-12
10% Tween-20	21-1038-12		

TECHNICAL INFORMATION

Storage	OPEN KIT IMMEDIATELY and store components at room temp, 4°C, -20°C, and -80°C as indicated (see User Manual corresponding to Kit Version 1). Stable for six months upon date of receipt.
Instructions for Use	See User Manual corresponding to Kit Version 1. Pair with CUTANA CDI Primers (14-1190-24rxn) for a full cells-to-libraries workflow.

REFERENCES

- [1] Buenrostro et al. *Nature Methods* (2013). PMID: 24097267
[2] Corces et al. *Nature Methods* (2017). PMID: 28846090
[3] Grandi et al. *Nature Protocols* (2022). PMID: 35478247

SAFETY INFORMATION

Item	Cat No	Chemical	Signal Word	Pictogram
Lysis Buffer	21-1039-12	NP-40	Warning	
2X Tagmentation Buffer	21-1040-12	DMF	Danger	  
Non-Hot Start 2X PCR Master Mix	15-1018-12	DMSO	Warning	
5% Digitonin	21-1004-12	Digitonin, DMSO	Danger	

CERTIFICATE OF ANALYSIS

Quality Control Metrics	Result
Functional Testing (ATAC-seq) Each lot of CUTANA™ ATAC-seq Kit is functionally validated by comparison to a previous lot and to established EpiCypher internal controls. ATAC-seq is performed on 50k native K562 nuclei using the CUTANA™ ATAC-seq Kit. Libraries are sequenced on an Illumina NextSeq2000 with paired-end sequencing (2x50 bp). Assessment is based on metrics including library fragment size distribution and yields, percent unique reads, percent mitochondrial reads, percent E. coli reads, fraction of reads in peaks (FRiP), insert size distribution, and TSS enrichment score.	Pass

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.