



A Hologic Company

D-Plex Single Indexes Module

Indexes modules for D-Plex Small RNA-seq Kit

Cat. No. C05030010 (Set A: 24 indexes, 24 rxns)
C05030011 (Set B: 24 indexes, 24 rxns)

USER GUIDE

Version 2 03_2021

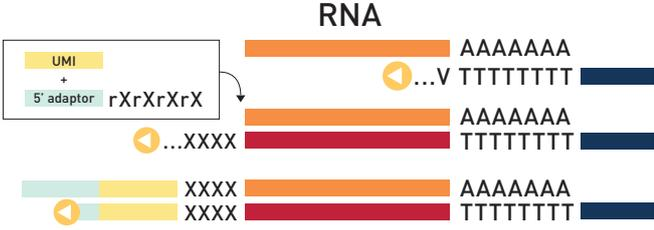
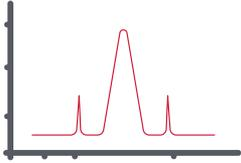
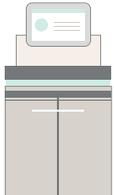


Please read D-Plex Small RNA-seq manual carefully before starting your experiment

Summary

Kit Method Overview	4
Introduction	5
Materials	7
Multiplexing Advices	10
Related Products	11

Kit Method Overview

			Time	Pause
STEP 1		RNA 	RNA sample preparation	5 min 
STEP 2		RNA 	RNA tailing	45 min
STEP 3		RNA 	Reverse transcription with template switching	>2h 
STEP 4		DNA 	PCR amplification	45 min 
STEP 5			Library purification	
STEP 6			Library clean-up and shaping	
STEP 7			Library quality control and quantification	

Introduction

The Diagenode D-Plex Small RNA-seq Library Preparation Kit is a tool designed for the study of the **small non-coding transcriptome**. The present kit incorporates the unique **D-Plex technology** to generate small RNA libraries for **Illumina sequencing**.

The D-Plex technology utilizes the innovative **capture and amplification by tailing and switching**, a ligation-free method for library preparation and offers key advantages such as:

- **Ultra-low input** capability of the library preparation
- Ease of use in a **one day, one tube protocol**
- **Higher library complexity** than most of other available library preparation kits for small RNA-sequencing

The library preparation protocol works on either total intact RNAs (RIN \geq 8) extracted and purified from a given sample or a small RNA fraction (<200nt), that might very well represent the circulating content of a **liquid biopsy-type of sample** (blood serum and plasma). The input requirements of the method are flexible and allow the user to perform the method within a wide range of RNA quantities going **from 10 pg** of small RNAs (<200nt) or circulating RNAs **up to 100 ng** of total RNAs.

The core of the technology relies on **ligation-free reactions** to attach the Illumina adaptors to both ends of the library construct. Therefore, the results generated with the D-Plex Small RNA-seq kit will vastly differ from those produced with ligase-based approach. For instance, the results generated with the D-Plex kit will encompass a **vast spectrum of small non-coding RNAs** (miRNAs, snoRNAs, snRNAs, piRNAs) whereas a ligase-based approach will enrich the sequencing library in 5'-P – 3'-OH RNAs, mainly mature miRNAs.

Diagenode therefore recommends having a **clear understanding of the scientific question** being asked in a given experiment before proceeding to a small RNA-seq library preparation as the choice of technology will strongly impact the end result.

For optimal workflow flexibility, the library preparation is available in both **unique dual index (UDI)** and **single index (SI)** configurations. The D-Plex Unique Dual Indexes Modules (C05030021 and C05030022) and the D-Plex Single Indexes Modules (C05030010 and C05030011) are available separately from the library preparation kit, providing PCR primers for library multiplexing up to 48. The use of UDI is highly recommended to mitigate errors introduced by read misassignment, including index hopping frequently observed with patterned flow cells such as Illumina's NovaSeq system.

An important addition to the D-Plex set of features is the use of **unique molecular identifiers (UMI)** to each transcript incorporated in the library. Given this new addition, it is now possible to exclude PCR duplicates from a set of reads, thus improving the transcript expression quantification.

Materials

Table 1: D-Plex Small RNA-seq Single Indexes Sequence - Set A (1-24)

D-Plex SI Reverse Primer Index #	PCR reverse primer sequence	Index
1	CAAGCAGAAGACGGCATAACGAGATCGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	ATCACG
2	CAAGCAGAAGACGGCATAACGAGATACATCGGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	CGATGT
3	CAAGCAGAAGACGGCATAACGAGATGCCTAAGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	TTAGGC
4	CAAGCAGAAGACGGCATAACGAGATTGGTCAAGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	TGACCA
5	CAAGCAGAAGACGGCATAACGAGATCACTGTGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	ACAGTG
6	CAAGCAGAAGACGGCATAACGAGATATTGGCGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	GCCAAT
7	CAAGCAGAAGACGGCATAACGAGATGATCTGGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	CAGATC
8	CAAGCAGAAGACGGCATAACGAGATTCAAGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	ACTTGA
9	CAAGCAGAAGACGGCATAACGAGATCTGATCGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	GATCAG
10	CAAGCAGAAGACGGCATAACGAGATAAGCTAGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	TAGCTT
11	CAAGCAGAAGACGGCATAACGAGATGTAGCCGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	GGCTAC
12	CAAGCAGAAGACGGCATAACGAGATTACAAGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	CTTGTA
13	CAAGCAGAAGACGGCATAACGAGATTTGACTGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	AGTCAA
14	CAAGCAGAAGACGGCATAACGAGATGGAAGTGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	AGTTCC
15	CAAGCAGAAGACGGCATAACGAGATTGACATGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	ATGTCA
16	CAAGCAGAAGACGGCATAACGAGATGGACGGGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	CCGTCC
17	CAAGCAGAAGACGGCATAACGAGATCTCTACGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	GTAGAG
18	CAAGCAGAAGACGGCATAACGAGATGCGGACGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	GTCCGC
19	CAAGCAGAAGACGGCATAACGAGATTTTACGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	GTGAAA
20	CAAGCAGAAGACGGCATAACGAGATGGCCACGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	GTGGCC
21	CAAGCAGAAGACGGCATAACGAGATCGAAACGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	GTTTCG
22	CAAGCAGAAGACGGCATAACGAGATCGTACGGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	CGTACG
23	CAAGCAGAAGACGGCATAACGAGATCCACTCGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	GAGTGG
24	CAAGCAGAAGACGGCATAACGAGATGCTACCGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	GGTAGC

(*) = phosphorothioate bond

Table 2: D-Plex Small RNA-seq Single Indexes Sequence - Set B (25-48)

D-Plex SI Reverse Primer Index #	PCR reverse primer sequence	Index
25	CAAGCAGAAGACGGCATAACGAGATATCAGTGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	ACTGAT
26	CAAGCAGAAGACGGCATAACGAGATGCTCATGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	ATGAGC
27	CAAGCAGAAGACGGCATAACGAGATAGGAATGTGACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	ATTCTT
28	CAAGCAGAAGACGGCATAACGAGATCTTTTGGTACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	CAAAAG
29	CAAGCAGAAGACGGCATAACGAGATTAGTTGGTACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	CAACTA
30	CAAGCAGAAGACGGCATAACGAGATCCGGTGGTACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	CACCGG
31	CAAGCAGAAGACGGCATAACGAGATATCGTGGTACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	CACGAT
32	CAAGCAGAAGACGGCATAACGAGATTGAGTGGTACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	CACTCA
33	CAAGCAGAAGACGGCATAACGAGATCGCCTGGTACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	CAGGCG
34	CAAGCAGAAGACGGCATAACGAGATGCCATGGTACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	CATGGC
35	CAAGCAGAAGACGGCATAACGAGATAAAATGGTACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	CATTTT
36	CAAGCAGAAGACGGCATAACGAGATTGTTGGGTACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	CCAACA
37	CAAGCAGAAGACGGCATAACGAGATATTCCGGTACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	CGGAAT
38	CAAGCAGAAGACGGCATAACGAGATAGCTAGGTACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	CTAGCT
39	CAAGCAGAAGACGGCATAACGAGATGTATAGGTACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	CTATAC
40	CAAGCAGAAGACGGCATAACGAGATTCTGAGGTACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	CTCAGA
41	CAAGCAGAAGACGGCATAACGAGATGTCGTCGTACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	GACGAC
42	CAAGCAGAAGACGGCATAACGAGATCGATTAGTACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	TAATCG
43	CAAGCAGAAGACGGCATAACGAGATGCTGTAGTACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	TACAGC
44	CAAGCAGAAGACGGCATAACGAGATATTATAGTACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	TATAAT
45	CAAGCAGAAGACGGCATAACGAGATGAATGAGTACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	TCATTC
46	CAAGCAGAAGACGGCATAACGAGATTCGGGAGTACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	TCCCGA
47	CAAGCAGAAGACGGCATAACGAGATCTTCGAGTACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	TCGAAG
48	CAAGCAGAAGACGGCATAACGAGATTGCCGAGTACTGGAGTTCAGACGTGTGCTCTTCCGATC*T	TCGGCA

(*) = phosphorothioate bond

Table 3: Module content

Component	Cap color	Quantity	Storage
D-Plex Forward Primer FP (x1)	Blue	240 µl	-20°C/-4°F
D-Plex Reverse Primer Index RP (x24)	Blue	30 µl each	-20°C/-4°F
RT Primer H (RTPH)	Purple	24 µl	-20°C/-4°F
RT Primer M (RTPM)	Purple	24 µl	-20°C/-4°F
Small Template Switching Oligo (STSO)	Purple	48 µl	-20°C/-4°F

Important Notice

The RT Primer H, RT Primer M and Small Template Switching Oligo components included in the D-Plex Small RNA-seq kit (C05030001) are only suitable for unique dual index (UDI) library construction.

For single index (SI) library construction, the RT Primer H, RT Primer M and Small Template Switching Oligo components suitable for SI library constructions are included in the D-Plex Single Indexes modules (C05030010 and C05030011). You should use the components corresponding to the desired – UDI or SI – library construction.

Multiplexing Advices

The D-Plex PCR reverse primers in Table 1 and 2 bear the TruSeq (Illumina) Small RNA adapters that can be used for library **multiplexing up to 48**.

In case of a multiplexing scenario, we recommend to follow Illumina's library pooling guidelines that are explained in Table 4 and submit the D-Plex libraries as TruSeq small RNA libraries to your sequencing provider.

Table 4: Multiplexing recommendations for the D-Plex Small RNA-seq SI library construction

Level of multiplexing	Option #	For index set A (#1-24)
2	1	RP#6 + RP#12
3	1	RP#1 + RP#3 + RP#7
	2	RP#2 + RP#4 + RP#8
	3	RP#16 + RP#17 + RP#18
	4	RP#13 + RP#17 + RP#23
	5	2-plex with any other index
4	1	RP#2 + RP#9 + RP#10 + RP#11
	2	RP#4 + RP#5 + RP#6 + RP#7
	3	3-plex with any other index
Level of multiplexing	Option #	For index set B (#25-48)
2	1	RP#37 + RP#45
3	1	RP#38 + RP#44 + RP#46
	2	RP#40 + RP#47 + RP#48
	3	2-plex with any other index
4	1	RP#37 + RP#39 + RP#42 + RP#43
	2	RP#37 + RP#38 + RP#45 + RP#46
	3	3-plex with any other index

Related Products

Product	Reference
D-Plex Small RNA-seq Kit	C05030001
D-Plex Unique Dual Indexes Module – Set A	C05030021
D-Plex Unique Dual Indexes Module – Set B	C05030022

FOR RESEARCH USE ONLY.

Not intended for any animal or human therapeutic or diagnostic use.

© 2021 Diagenode SA. All rights reserved. No part of this publication may be reproduced, transmitted, transcribed, stored in retrieval systems, or translated into any language or computer language, in any form or by any means: electronic, mechanical, magnetic, optical, chemical, manual, or otherwise, without prior written permission from Diagenode SA (hereinafter, "Diagenode"). The information in this guide is subject to change without notice. Diagenode and/or its affiliates reserve the right to change products and services at any time to incorporate the latest technological developments. Although this guide has been prepared with every precaution to ensure accuracy, Diagenode and/or its affiliates assume no liability for any errors or omissions, nor for any damages resulting from the application or use of this information. Diagenode welcomes customer input on corrections and suggestions for improvement.

NOTICE TO PURCHASER LIMITED LICENSE

The information provided herein is owned by Diagenode and/or its affiliates. Subject to the terms and conditions that govern your use of such products and information, Diagenode and/or its affiliates grant you a nonexclusive, nontransferable, non-sublicensable license to use such products and information only in accordance with the manuals and written instructions provided by Diagenode and/or its affiliates. You understand and agree that except as expressly set forth in the terms and conditions governing your use of such products, that no right or license to any patent or other intellectual property owned or licensable by Diagenode and/or its affiliates is conveyed or implied by providing these products. In particular, no right or license is conveyed or implied to use these products in combination with any product not provided or licensed to you by Diagenode and/or its affiliates for such use. Limited Use Label License: Research Use Only The purchase of this product conveys to the purchaser the limited, non-transferable right to use the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact info@diagenode.com.

TRADEMARKS

The trademarks mentioned herein are the property of Diagenode or their respective owners. Bioanalyzer is a trademark of Agilent Technologies, Inc. Agencourt and AMPure® are registered trademarks of Beckman Coulter, Inc. Illumina® is a registered trademark of Illumina® Inc; Qubit is a registered trademark of Life Technologies Corporation.

www.diagenode.com