

D-Plex Unique Dual Indexes

For use with the D-Plex small RNA-seq kit

Set A: 24 UDIs, 24 rxns

Cat. No. C05030021

Set B: 24 UDIs, 24 rxns

Cat. No. C05030022

Set C: 24 UDIs, 24 rxns

Cat. No. C05030023

Set D: 24 UDIs, 24 rxns

Cat. No. C05030024

Content

Introduction	4
Materials	5
Multiplexing Advices.....	9
Related Products.....	10
Revision History.....	10



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

Introduction

The Hologic Diagenode D-Plex Unique Dual Indexes are an essential component of the D-Plex RNA sequencing library preparation workflow. These unique dual indexes (UDI) are designed and validated for use with the D-Plex Small RNA-seq Kit for Illumina sequencing platform (Cat. No. C05030001).

The D-Plex technology utilizes the innovative **Capture and Amplification by Tailing and Switching (CATS)**, a ligation-free method for library preparation, and offers key advantages such as:

- **Low to ultra-low input** capability
- Ease of use in a **one-day, one-tube protocol**
- **High performance** and **library complexity**

Each D-Plex Unique Dual Indexes module – Set A, B, C, or D – includes 24 primer pairs with **unique dual barcodes** (unique i5 and i7 indexes) for library multiplexing of up to 24. Four sets – Set A, B, C, and D – can be used simultaneously for library **multiplexing of up to 96**. The use of UDIs 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.

Materials

Table 1. D-Plex Unique Dual Indexes Sequence – Set A (1-24)

D-Plex Primer UDI #	PCR Reverse Primer		PCR Forward Primer	
	i7 Bases for sample sheet	i5 bases for sample sheet Forward Strand Workflow NovaSeq 6000 v1.0, MiSeq, HiSeq 2000/2500	i5 bases for sample sheet Reverse Strand Workflow NovaSeq 6000 v1.5, iSeq, MiniSeq, NextSeq, HiSeq 3000/4000	
1	CCGCGGTT	AGCGCTAG	CTAGCGCT	
2	TTATAACC	GATATCGA	TCGATATC	
3	GGA CTTGG	CGCAGACG	CGTCTGCG	
4	AAGTCCAA	TATGAGTA	TACTCATA	
5	ATCCACTG	AGGTGCGT	ACGCACCT	
6	GCTTGTC A	GAACATAC	GTATGTTC	
7	CAAGCTAG	ACATAGCG	CGCTATGT	
8	TGGATCGA	GTGCGATA	TATCGCAC	
9	AGTTCAGG	CCAACAGA	TCTGTTGG	
10	GACCTGAA	TTGGTGAG	CTCACCAA	
11	TCTCTACT	CGCGGTTC	GAACCGCG	
12	CTCTCGTC	TATAACCT	AGGTTATA	
13	CCAAGTCT	AAGGATGA	TCATCCTT	
14	TTGGACTC	GGAAGCAG	CTGCTTCC	
15	GGCTTAAG	TCGTGACC	GGTCACGA	
16	AATCCGGA	CTACAGTT	AACTGTAG	
17	TAATACAG	ATATTCAC	GTGAATAT	
18	CGGCGTGA	GCGCCTGT	ACAGGCGC	
19	ATGTAAGT	ACTCTATG	CATAGAGT	
20	GCACGGAC	GTCTCGCA	TGCGAGAC	
21	GGTACCTT	AAGACGTC	GACGTCTT	
22	AACGTTCC	GGAGTACT	AGTACTCC	
23	GCAGAATT	ACCGGCCA	TGGCCGGT	
24	ATGAGGCC	GTTAATTG	CAATTAAC	

Table 2. D-Plex Unique Dual Indexes Sequence – Set B (25-48)

D-Plex Primer UDI #	PCR Reverse Primer	PCR Forward Primer	
	i7 Bases for sample sheet	i5 bases for sample sheet Forward Strand Workflow NovaSeq 6000 v1.0, MiSeq, HiSeq 2000/2500	i5 bases for sample sheet Reverse Strand Workflow NovaSeq 6000 v1.5, iSeq, MiniSeq, NextSeq, HiSeq 3000/4000
25	ACTAAGAT	AACCGCGG	CCGCGGTT
26	GTCGGAGC	GGTTATAA	TTATAACC
27	CTTGGTAT	CCAAGTCC	GGACTTGG
28	TCCAACGC	TTGGACTT	AAGTCCAA
29	CCGTGAAG	CAGTGGAT	ATCCACTG
30	TTACAGGT	TGACAAGC	GCTTGTC A
31	GGCATTCT	CTAGCTTG	CAAGCTAG
32	AATGCCTC	TCGATCCA	TGGATCGA
33	TACCGAGG	CCTGAACT	AGTTCAGG
34	CGTTAGAA	TTCAGGTC	GACCTGAA
35	AGCCTCAT	AGTAGAGA	TCTCTACT
36	GATTCTGC	GACGAGAG	CTCTCGTC
37	TCGTAGTG	AGACTTGG	CCAAGTCT
38	CTACGACA	GAGTCCAA	TTGGACTC
39	TAAGTGGT	CTTAAGCC	GGCTTAAG
40	CGGACAAC	TCCGGATT	AATCCGGA
41	ATATGGAT	CTGTATTA	TAATACAG
42	GCGCAAGC	TCACGCCG	CGGCGTGA
43	AAGATACT	ACTTACAT	ATGTAAGT
44	GGAGCGTC	GTCCGTGC	GCACGGAC
45	ATGGCATG	AAGGTACC	GGTACCTT
46	GCAATGCA	GGAACGTT	AACGTTCC
47	GTTCCAAT	AATTCTGC	GCAGAATT
48	ACCTTGGC	GGCCTCAT	ATGAGGCC

Table 3. D-Plex Unique Dual Indexes Sequence – Set C (49-72)

D-Plex Primer UDI #	PCR Reverse Primer	PCR Forward Primer	
	i7 Bases for sample sheet	i5 bases for sample sheet Forward Strand Workflow NovaSeq 6000 v1.0, MiSeq, HiSeq 2000/2500	i5 bases for sample sheet Reverse Strand Workflow NovaSeq 6000 v1.5, iSeq, MiniSeq, NextSeq, HiSeq 3000/4000
49	ATATCTCG	ATCTTAGT	ACTAAGAT
50	GCGCTCTA	GCTCCGAC	GTCGGAGC
51	AACAGGTT	ATACCAAG	CTTGGTAT
52	GGTGAACC	GCGTTGGA	TCCAACGC
53	CAACAATG	CTTCACGG	CCGTGAAG
54	TGGTGGCA	TCCTGTAA	TTACAGGA
55	AGGCAGAG	AGAATGCC	GGCATTCT
56	GAATGAGA	GAGGCATT	AATGCCTC
57	TGCGGCGT	CCTCGGTA	TACCGAGG
58	CATAATAC	TTCTAACG	CGTTAGAA
59	GATCTATC	ATGAGGCT	AGCCTCAT
60	AGCTCGCT	GCAGAATC	GATTCTGC
61	CGGAACTG	CACTACGA	TCGTAGTG
62	TAAGGTCA	TGTCGTAG	CTACGACA
63	TTGCCTAG	ACCACTTA	TAAGTGGT
64	CCATTCGA	GTTGTCCG	CGGACAAC
65	ACACTAAG	ATCCATAT	ATATGGAT
66	GTGTCGGA	GCTTGCGC	GCGCAAGC
67	TTCCTGTT	AGTATCTT	AAGATACT
68	CCTTCACC	GACGCTCC	GGAGCGTC
69	GCCACAGG	CATGCCAT	ATGGCATG
70	ATTGTGAA	TGCATTGC	GCAATGCA
71	ACTCGTGT	ATTGGAAC	GTTCCAAT
72	GTCTACAC	GCCAAGGT	ACCTTGGC

Table 4. D-Plex Unique Dual Indexes Sequence – Set D (73-96)

D-Plex Primer UDI #	PCR Reverse Primer	PCR Forward Primer	
	i7 Bases for sample sheet	i5 bases for sample sheet Forward Strand Workflow NovaSeq 6000 v1.0, MiSeq, HiSeq 2000/2500	i5 bases for sample sheet Reverse Strand Workflow NovaSeq 6000 v1.5, iSeq, MiniSeq, NextSeq, HiSeq 3000/4000
73	CAATTAAC	CGAGATAT	ATATCTCG
74	TGGCCGGT	TAGAGCGC	GCGCTCTA
75	AGTACTCC	AACCTGTT	AACAGGTT
76	GACGTCTT	GGTTCACC	GGTGAACC
77	TGCGAGAC	CATTGTTG	CAACAATG
78	CATAGAGT	TGCCACCA	TGGTGGCA
79	ACAGGCGC	CTCTGCCT	AGGCAGAG
80	GTGAATAT	TCTCATTC	GAATGAGA
81	AACTGTAG	ACGCCGCA	TGCGGCGT
82	GGTCACGA	GTATTATG	CATAATAC
83	CTGCTTCC	GATAGATC	GATCTATC
84	TCATCCTT	AGCGAGCT	AGCTCGCT
85	AGGTTATA	CAGTTCCG	CGGAACTG
86	GAACCGCG	TGACCTTA	TAAGGTCA
87	CTCACCAA	CTAGGCAA	TTGCCTAG
88	TCTGTTGG	TCGAATGG	CCATTCGA
89	TATCGCAC	CTTAGTGT	ACACTAAG
90	CGCTATGT	TCCGACAC	GTGTGCGA
91	GTATGTTC	AACAGGAA	TTCCTGTT
92	ACGCACCT	GGTGAAGG	CCTTCACC
93	TACTCATA	CCTGTGGC	GCCACAGG
94	CGTCTGCG	TTCACAAT	ATTGTGAA
95	TCGATATC	ACACGAGT	ACTCGTGT
96	CTAGCGCT	GTGTAGAC	GTCTACAC

Table 5. Module Content

Component	Cap color	Quantity	Storage temperature
D-Plex Primer UDI (x24)	Black	20 µl each (= 1 rxn/UDI)	-20°C/-4°F

Multiplexing Advices

The D-Plex PCR primers in Tables 1 to 4 bear TruSeq (Illumina) HT adapters, which can be used for library **multiplexing of up to 96**.

In a multiplexing scenario, we recommend following the numerical order of the indexes and submitting the D-Plex libraries as TruSeq HT libraries to your sequencing provider. For example, if 8 samples need to be multiplexed together, one can simply select a subset of 8 indexes following the consecutive order from the D-Plex Unique Dual Indexes module – Set A or from the D-Plex Unique Dual Indexes module – Set B, C, or D.

Related Products

Product	Cat. No.
D-Plex Small RNA-seq Kit	C05030001

Revision history

Version	Date of modification	Description of modifications
Version 2 04_2026	April 2026	Removal of obsolete references. Correction of a typo.
Version 2 09_2022	September 2022	Addition of the information related to the Sets C & D
Version 1 03_2021	March 2021	Manual Creation

FOR RESEARCH USE ONLY.

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

© 2026 Hologic Diagenode S.A. 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 Hologic Diagenode. The information in this guide is subject to change without notice. Hologic 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, Hologic 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. Hologic Diagenode welcomes customer input on corrections and suggestions for improvement.

NOTICE TO PURCHASER LIMITED LICENSE

The information provided herein is owned by Hologic Diagenode and/or its affiliates. Subject to the terms and conditions that govern your use of such products and information, Hologic 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 Hologic 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 Hologic 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 Hologic 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 Hologic 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.

diagenode.com